

# Attune™ Cytometric Software

## USER GUIDE

For data acquisition and analysis using the Attune™ NxT and  
Attune™ CytPix™ Flow Cytometers

Publication Number MAN0026553

Revision D



Revision history: MAN0026553 D (English)

Revision	Date	Description
D	14 October 2024	Add information about Baseline Functional Response, add information about local user access when SAE is enabled, update all relevant screenshots.
C.0	30 January 2024	Update the Off-plot gates section and add info about off-plot gate indicators; add Maintenance tab to System Log dialog, update Instrument and Workspace Ribbon tabs; update Experiment Explorer context menus to add Remove Image Processing Data; update File Dialogs to add Compression Level options; update Appendix D: SAE Administrator Console - Roles, System, and Audit History tabs.
B.0	27 April 2023	Add Process Images dialog and Attune Image Processing Dashboard chapters; add Attune Cytometric SW Image processing parameters appendix; update Ribbon tabs, Image View, Experiment Explorer, Customize panel, Image Capture Settings panel, Options dialog, File dialogs, Dialogs, Attune Database Utility chapters and the Data Management appendix with information about image processing and data processing; convert the UG to CCMS and change the formatting and style.
A.0	25 April 2022	Change Pub. No. from 100024236 to MAN0026553, add Autogating section to WS tab and to Gate options, add Autogate controls to Customize Gates panel, update UI screens to include Autogating tools and updated icons, update Cloud sign in and Enable SAE mode workflows.

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

**TRADEMARKS:** All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2015-2024 Thermo Fisher Scientific Inc. All rights reserved.



# Contents

■ About this guide .....	28
Conventions .....	28
Text and keyboard conventions .....	28
Clicking .....	28
User attention symbols .....	29
Other Attune™ user guides .....	29
Safety information .....	30
Safety alert words .....	30
SDSs .....	30
■ CHAPTER 1 Attune™ Cytometric Software .....	31
Attune™ Cytometric Software overview .....	31
About the software .....	31
System requirements .....	32
Supported files .....	32
Software layout and ribbon bar menu system .....	32
Software layout .....	32
Ribbon bar menu system .....	32
Context menus .....	33
Keyboard shortcuts .....	33
Attune™ Cytometric Software installation .....	33
Guidelines for installation .....	34
Install the Attune™ Cytometric Software .....	34
Update the instrument firmware .....	35
■ CHAPTER 2 Startup, Login, and Main Menu .....	36
Main application startup .....	36
Main application startup .....	36
End user license agreement (EULA) .....	36
SAE module .....	37
Login screen .....	37
Sign in – standard mode .....	38
Sign in – SAE mode .....	38
SAE account rules .....	40

Invalid username or password .....	40
Expired password .....	41
Account types .....	42
Default local Attune™ accounts .....	42
Main Menu .....	46
Log Out .....	47
Main Menu ribbon bar .....	47
Experiment Explorer .....	48
<b>■ CHAPTER 3 Attune™ Desktop .....</b>	<b>49</b>
Attune™ Desktop overview .....	49
Attune™ Desktop layout .....	50
Application title bar .....	51
<b>Quick Access toolbar</b> .....	52
Default buttons .....	52
Sign in to Cloud .....	53
Ribbon bar .....	55
Ribbon tabs .....	55
Ribbon context menu .....	55
<b>Main application workspace</b> .....	56
Application area .....	56
Application panels .....	57
Docking locations .....	58
Docking and undocking panels .....	59
Docking and undocking tabs .....	60
Floating panels .....	60
Auto Hide .....	60
<b>Application status bar</b> .....	63
Instrument status and alerts .....	63
Camera status .....	66
Connect status .....	66
Tube and Plate modes .....	66
Time .....	68
Coordinates .....	68
Workspace and Instrument Settings .....	69
Acquisition status .....	69
Analysis status .....	69
Size slider .....	70

■ CHAPTER 4 Ribbon tabs .....	71
Overview .....	71
Ribbon tabs .....	71
File tab .....	72
File menu .....	72
Home tab .....	74
Create group .....	74
Templates group .....	74
General group .....	75
Clipboard group .....	75
Editing group .....	76
Refresh group .....	76
View tab .....	77
Experiment group .....	77
Show Panels group .....	78
Show Colors .....	79
Other Options group .....	80
Workspace tab .....	81
Save as Default Workspace .....	81
Plots group .....	82
Gating Tools group .....	83
Autogating Tools group .....	84
Cell Image Analysis group .....	85
Other group .....	85
Size and Position group .....	86
Instrument tab .....	88
Setup group .....	88
System History group .....	89
Functions group .....	90
Service group .....	93
Automation group .....	94
Experiment tab .....	95
Experiment limits .....	95
Experiment tab controls .....	95
Hot keys .....	98
Compensation tab .....	99
Setup group .....	99
Apply group .....	100
Adjustment group .....	101
Statistics tab .....	103
Statistics tab behavior .....	103
Results group .....	105
Tools group .....	106
General group .....	107

Event Statistics group .....	108
Intensity group .....	108
Variation group .....	109
Overlay tab .....	110
Create group .....	110
Gating Tools group .....	111
Galleries View group .....	111
Image Settings tab .....	112
Setup group .....	112
Analysis group .....	113
Image Processing group .....	113
SAE tab .....	114
Signing group .....	114
Other group .....	115
■ <b>CHAPTER 5</b> Workspace view .....	<b>116</b>
Overview .....	116
Workspace behavior .....	117
Workspace selection .....	118
Workspace levels .....	118
Workspace view label .....	118
Workspace selection menu .....	118
Workspace objects .....	120
Create Workspace objects .....	120
Select Workspace objects .....	121
Resize Workspace objects .....	123
Move Workspace objects .....	124
Delete, format, or duplicate Workspace objects .....	124
Paste Workspace objects .....	125
Workspace view layout .....	126
Resize the Workspace view .....	126
Workspace view modes .....	126
Freeform mode .....	126
Auto Layout mode .....	127
Previews .....	132
Overview .....	132
Access the Preview panel .....	132
Preview panel title bar .....	133
Preview plots .....	134
Resize Preview plots .....	134
Add Plots to Workspace – Single Plot selection .....	134
Add Plots to Workspace – Group selection .....	135
Mouse over behavior .....	136

Plots .....	138
Overview .....	138
Histogram plot .....	139
Dot plot .....	140
Density plot .....	141
Precedence density plot .....	142
Adjust plots .....	143
Gates .....	145
Overview .....	145
Create and delete gates .....	145
Gate name .....	146
Gate color .....	147
Move and resize gates .....	148
Rectangular Gate .....	149
Oval Gate .....	150
Polygon Gate .....	151
Quadrant Gate .....	152
Bent Quadrant Gate .....	153
Histogram Gate .....	155
Contour Autogate .....	156
Ellipse autogate .....	159
Autogate gating rules .....	162
Derived gate .....	163
Off-plot gates .....	164
Gate behavior .....	173
Workspace statistics .....	175
Overview .....	175
Workspace statistics .....	176
Plot statistics (single plot) .....	177
Plot statistics (multiple plots) .....	177
Workspace images and text .....	178
Workspace images .....	178
Workspace text .....	178
Image backgating .....	179
Cell image container .....	181
Overview .....	181
Population menu .....	182
Navigation .....	183
Measure Image tool .....	184
Workspace view context menu .....	185
Overview .....	185
Insert .....	185
Paste and Select all .....	186
Send All Plots To Overlay .....	187
Plot title hyperlink .....	188

Plot context menu .....	189
Overview .....	189
Customize .....	189
Scale .....	190
Set Population .....	191
Add gate .....	192
Insert plot statistics .....	192
Swap Axis .....	192
Cut, Copy, and Delete .....	192
Send to overlay .....	193
Save as .....	193
Gate context menu .....	194
Overview .....	194
Customize .....	194
Create daughter plot .....	195
Export FCS file .....	195
Cut, Copy, and Delete .....	196
Edit gates .....	196
Plot axis context menu .....	197
Overview .....	197
Customize .....	197
Scale .....	198
Parameter .....	199
Set Population .....	201
Add gate .....	202
Insert plot statistics .....	202
Swap Axis .....	202
Cut, Copy, Paste, and Delete .....	202
Send to overlay .....	203
Save as .....	204
Plot parameter hyperlink .....	205
Statistics context menu .....	207
Overview .....	207
Customize .....	207
Cut, Copy, and Delete .....	207
Export stats .....	207
Image context menu .....	208
Overview .....	208
Cut, Copy, and Delete .....	208
Text box context menu .....	208
Overview .....	208
Customize .....	209
Cut, Copy, and Delete .....	209

Text box text context menu .....	209
Overview .....	209
Cut, Copy, and Delete .....	209
Cell Image Container context menu .....	210
Overview .....	210
Customize .....	210
Cut, Copy, and Delete .....	210
Measure .....	210
<b>■ CHAPTER 6 Heat Map View .....</b>	<b>211</b>
Overview .....	211
Display of samples .....	212
Overview .....	212
Tube Experiment .....	213
Plate Experiment .....	214
Navigation and selection .....	215
Set up an Experiment .....	216
Experiment/Group labels .....	216
Compensation wells .....	216
Manual wells .....	217
Run Protocol .....	217
Legend .....	218
Analyze results .....	220
Display of results .....	220
Tooltips .....	220
Indicators .....	221
<b>■ CHAPTER 7 Sample List view .....</b>	<b>222</b>
Overview .....	222
Sample List properties .....	224
Edit values in the Sample List .....	224
Rules for editing the Sample List .....	224
Navigation in the Sample List with the keyboard .....	225
Selection in the Sample List .....	225
Multiselection in the Sample List .....	226
Copy and Paste .....	226
Print the Sample List .....	227
Sample List context menu .....	227
Sample List context menus .....	227

■	<b>CHAPTER 8 Results view</b>	<b>228</b>
	Overview	228
	Results table	229
	Results table	229
	Primary table	229
	Select rows	230
	Group results	230
	Group results by column headers	230
	Add and remove column headers	231
	Moving versus copying headers	232
	Examples of grouped results	232
	Results table column header context menu	233
■	<b>CHAPTER 9 Overlays</b>	<b>234</b>
	Overview	234
	Overlays tab	234
	Overlays area	235
	Galleries area	236
	Overlay plots	237
	Overlay plot modes	237
	Overlay plot characteristics	238
	Gallery plots	241
	Gallery plot characteristics	241
	Control plot	241
	Overlay calculations	243
	Display of statistics	244
	Gates	245
	Overlay properties	246
	Overlay settings	246
	Add plots to Overlays	247
	Compare Plots with different Scale and Range settings	247
	Select plots	248
	Move and copy plots between Overlays	250
	Copy plot images	251
	Resize plots	251
	Overlay plot links to Source plot	251
	Auto-Refresh	251
	Overlay plot context menu	253
	Overview	253
	Overlay plot context menu options	253



Gallery plot context menu .....	256
Overview .....	256
Gallery plot context menu options .....	256
Overlay Builder .....	260
Overview .....	260
Select and order samples .....	262
Select Plots .....	264
Page Setup .....	266
Print overlays .....	268
<b>■ CHAPTER 10 Image View .....</b>	<b>269</b>
Overview .....	269
Image view tab .....	270
Active image .....	271
Image gallery .....	272
Navigation and selection .....	273
Image capture gate .....	276
Image backgating .....	278
Measure image tool .....	280
Image view context menus .....	282
Active image context menu .....	282
Image Gallery context menu .....	283
Measure image tool context menu .....	284
<b>■ CHAPTER 11 Experiment Explorer .....</b>	<b>285</b>
Overview .....	285
Experiment Explorer icons .....	286
Files view .....	287
Organization .....	287
Storage and Size .....	288
Experiment hierarchy .....	291
Plate hierarchy .....	291
Tube hierarchy .....	292
Experiment Explorer behavior .....	294
Selection indicators .....	294
Selection .....	294
Drag and drop .....	295
Instrument Settings (IS) badge identification .....	297
Drag-and-drop from Experiment Explorer to Overlays .....	298

Drag-and-drop FCS files .....	298
Display ID .....	299
Experiment Explorer context menus .....	301
Overview .....	301
Validation of name fields .....	301
User and Shared folder context menus .....	302
Plate context menu .....	304
Experiment context menu .....	310
Group context menu .....	318
Sample context menu .....	324
Compensation node context menu .....	331
Compensation control context menu .....	335
Deleted items dialog .....	339

■ **CHAPTER 12** Collection panel ..... 340

Overview .....	340
Acquisition status .....	344
Acquisition status indicators .....	344
Sample remaining indicator .....	345
Collection controls .....	345
Overview .....	345
Collection controls – Tube and Manual wells .....	346
Overview of collection controls for tube and manual well samples .....	346
Buttons displayed .....	347
Collection controls – Plate .....	350
Overview of plate collection controls .....	350
Buttons displayed .....	352
Button availability .....	356
Current well indicator .....	356
Set up compensation workflow .....	357
Collect and display options .....	358
Collect group overview .....	358
Collect options .....	359
Record and save dialogs .....	364
Run protocol .....	368
Overview .....	368
Settings buttons .....	371
Optimize for high throughput collection .....	372
Set Up Comp Options .....	373
Flow/Recording Flow Options .....	374
Stop Options .....	377
Other options .....	379
Rinse reminder dialog .....	381

■ CHAPTER 13 Instrument settings panel .....	382
Overview .....	382
General panel properties .....	383
Default and load options .....	383
Overview .....	383
Automatic update option .....	383
Set as experiment IS .....	384
Save as Default split button .....	384
Parameters .....	385
Overview .....	385
Configuration dropdown .....	387
Update all samples .....	388
Parameter selector .....	389
Target textbox .....	390
Label textbox .....	391
Notification badge .....	391
Threshold .....	392
Overview .....	392
Logic selector .....	393
Channel selector .....	393
Threshold adjustment slider bar .....	394
Threshold textbox .....	394
Exclude coincident events check box .....	395
Voltage .....	396
Overview .....	396
Voltage adjustment slider bar .....	397
Voltage value textbox .....	397
Advanced™ settings .....	398
Overview .....	398
Width threshold setting .....	399
Area scaling factor (ASF) .....	400
Window extension setting .....	401
Notification button .....	401
System instrument settings .....	401
Overview .....	401
Area scaling factor (ASF) .....	403
Window extension setting .....	404
Notification icon .....	404

- **CHAPTER 14 Heat map setup panel** ..... 405
  - Overview ..... 405
  - Apply, save, and load options ..... 407
    - Overview ..... 407
    - Apply changes ..... 407
    - Set as default split button ..... 408
  - Experiment ..... 408
    - Overview ..... 408
    - Experiment dropdown ..... 409
    - Experiment color button ..... 409
    - Statistic dropdown ..... 410
    - Gate dropdown ..... 411
    - Parameter dropdown ..... 411
  - Display mode ..... 412
    - Overview ..... 412
    - Display mode options ..... 412
  - Sample mode – threshold ..... 414
    - Overview ..... 414
    - Transition colors ..... 415
    - Transition values ..... 415
    - Create additional transition points ..... 416
  - Sample mode – heat map ..... 417
    - Overview ..... 417
    - Transition colors ..... 418
    - Transition values ..... 418
    - Create additional transition points ..... 419
- **CHAPTER 15 Customize panel** ..... 420
  - Overview ..... 420
    - Customize panel behavior ..... 420
  - Customize plot options ..... 421
    - Overview ..... 421
    - Plot types ..... 421
    - Customize panel behavior for plots ..... 424
      - General group ..... 427
      - General group – plot-specific options ..... 434
      - X-axis group ..... 436
      - Y-axis group ..... 441
      - Text group (plots only) ..... 442
  - Customize text box options ..... 444
    - Text box options ..... 444
  - Customize statistics box options ..... 445
    - Statistics box options ..... 445

Customize gate options .....	447
Overview .....	447
Customize panel behavior for gates .....	447
Gate type group .....	448
Gate type buttons .....	454
Name options .....	459
Coordinates group .....	461
Autogating options .....	464
Customize experiment (Plate or Tube) options .....	466
Overview .....	466
Customize panel for plate experiments .....	467
Customize panel behavior for tubes .....	468
Plate information .....	468
Experiment information .....	470
Group information .....	471
Sample information .....	473
Customize Overlay options .....	474
Overview .....	474
General options .....	475
3D Options .....	479
Calculations .....	481
Text .....	483
Customize Image options .....	484
Overview .....	484
Image Adjustment .....	485
Mask Settings .....	489
SAE object properties .....	490
Overview .....	490
<b>■ CHAPTER 16 Image Capture Settings panel .....</b>	<b>491</b>
Overview .....	491
Settings options .....	493
Automatically update .....	493
Set as Experiment Level .....	493
Set as Default .....	493
Capture Settings .....	494
Image Gate .....	494
Image Frequency .....	495
Total Images and Storage Size .....	495
Camera settings .....	497
Focus .....	497
Illumination .....	498

Image Size and Position Settings .....	499
Size and Position .....	500
Advanced .....	502
<b>■ CHAPTER 17 FCS information panel .....</b>	<b>503</b>
Overview .....	503
FCS file information .....	505
Sample information group .....	505
Parameters group .....	507
Compensation group .....	508
System information group .....	509
Data format group .....	510
Other group .....	511
<b>■ CHAPTER 18 Filter configuration .....</b>	<b>512</b>
Filter configuration (FC) module .....	512
Overview .....	512
Launch the FC module .....	513
Configuration ribbon tab .....	513
General group .....	514
Configuration management group .....	514
Edit filters dialog .....	519
View group .....	524
Default filter label list .....	525
Filter configuration view .....	526
Overview .....	526
Default instrument configurations .....	527
Configuration options .....	537
Overview .....	537
Hardware/Virtual laser configuration .....	538
<b>■ CHAPTER 19 Performance testing .....</b>	<b>539</b>
Overview .....	539
Baseline calculation and Performance testing .....	539
Performance tracking workflow (when to run) .....	539
Baseline calculation and Performance test data .....	540
Baseline Functional Response (BFR) .....	540
Baseline and Performance Test (BL/PT) module .....	541
Open the BL/PT module .....	541
Set up and run a Baseline .....	542
Baseline setup screen .....	542
Set up and run a baseline .....	544
Baseline screen .....	545

Completion of Baseline .....	547
Baseline completion without errors .....	547
Baseline completion with errors .....	547
Run a Performance test .....	548
Performance test setup screen .....	548
Set up and run a Performance test .....	550
Performance test screen .....	552
Completion of Performance test .....	553
Performance test completion without errors .....	553
Performance test completion with errors .....	554
Performance test reports .....	554
<b>■ CHAPTER 20 Performance Test reports .....</b>	<b>555</b>
Overview .....	555
Performance tests .....	555
Baseline calculation and Performance test data .....	556
Performance test reports .....	556
Performance History Report .....	557
View the Performance History Report .....	557
Results plot .....	559
Print the performance history report .....	560
Current PT results .....	561
View the Current PT Results .....	561
Current PT Result report .....	562
PT Results table .....	563
Print the PT results report .....	564
Levey-Jennings Report .....	565
View the Levey-Jennings Report .....	565
Levey-Jennings plots .....	566
Print the report .....	569
Baseline Report .....	572
View the Baseline Test Results .....	572
Baseline Results .....	573
Baseline Results table .....	574
Print the Baseline Results report .....	575
Baseline Functional Response (BFR) Report .....	576
View the Baseline Functional Response (BFR) Report .....	576
Print the Baseline Functional Response (BFR) report .....	578

- **CHAPTER 21 Compensation** ..... 579
  - Overview ..... 579
    - Compensation ..... 579
    - Compensation controls ..... 579
  - Compensation setup ..... 580
    - Initiate compensation setup ..... 580
    - Compensation setup dialog ..... 581
  - Compensation control source – Tubes ..... 581
    - Measurement ..... 582
    - Select background fluorescence mode ..... 583
    - Select compensation parameters ..... 584
    - OK and cancel buttons ..... 584
  - Compensation control source – wells ..... 585
    - Measurement ..... 586
    - Select background fluorescence mode ..... 586
    - Select compensation parameters ..... 588
    - OK and cancel buttons ..... 590
  - Compensation workspace ..... 590
    - Overview ..... 590
    - Customize the compensation workspace ..... 590
    - Plot context menu ..... 592
    - Plot axis context menu ..... 592
    - Gate context menu ..... 594
    - Compensation workspace for histogram view ..... 594
    - Gating plot ..... 596
    - Histogram compensation plot without negative gate ..... 597
    - Histogram compensation plot with negative gate ..... 598
    - Unstained control workspace ..... 599
  - Matrix dialog ..... 599
    - Overview ..... 599
    - Matrix dialog table ..... 600
    - OK, reset, and cancel buttons ..... 601
  - On-Plot compensation adjustment ..... 602
    - Overview ..... 602
    - Manual on-plot compensation adjustment ..... 603
  - Compensation acquisition workflow ..... 603
    - Overview ..... 603
    - Compensation acquisition from tubes ..... 603
    - Compensation acquisition from wells ..... 604
    - Compensation of instrument data ..... 604



■	<b>CHAPTER 22</b>	<b>New Experiment and Group dialogs</b>	<b>606</b>
		New Experiment dialog	606
		Open the New Experiment dialog	606
		Create a Tube-only experiment	607
		Create a plate experiment	609
		New experiment from template dialog	613
		Launch the new experiment from template dialog	614
		Select template	615
		Create experiment	618
		New group dialog	619
		Launch the new group dialog	619
		Create a group	620
■	<b>CHAPTER 23</b>	<b>Process Images dialog</b>	<b>622</b>
		Overview	622
		Open the Process Images dialog	623
		Select Samples	624
		Select Population	626
		Image Processing Options	628
		Image processing queue	630
		Completion of image processing	632
		Reprocess images	634
		Remove Image Processing Data dialog	635
■	<b>CHAPTER 24</b>	<b>Attune™ Image Processing Dashboard</b>	<b>636</b>
		Attune™ Image Processing Dashboard	636
		View image processing queue, status, and progress	638
		Reprioritize the image processing queue	638
		Cancel image processing requests	638
		Pause or resume image processing	638
■	<b>CHAPTER 25</b>	<b>Options dialogs</b>	<b>639</b>
		Overview	639
		General	641
		Overview	641
		New Experiment Defaults	642
		Naming Options	644
		Colors and Themes	647
		Overview	647
		Colors and Themes options	647

Fonts and Styles .....	649
Overview .....	649
Plot Title and Axis Label Style .....	650
Text Box Style .....	651
Previews .....	652
Plot Options .....	653
Overview .....	653
Ungated Color .....	653
Histogram Options .....	654
Dot Plot Options .....	654
Density Plot Options .....	655
Precedence Density Plot Options .....	655
Gate Options .....	656
Overview .....	656
Autogating .....	657
Gate Style .....	659
Gate Label Style .....	659
Gate Label Naming .....	660
Export Options .....	662
Overview .....	662
Plot Export Options .....	663
Plot export preview .....	663
Default Plot Image Format .....	664
Export Size (scaled to width) .....	664
Stats Options .....	665
Overview .....	665
Header .....	666
Statistics .....	666
Stats Box Style .....	667
Decimal Settings .....	667
Keyword Options .....	669
Overview .....	669
Keyword Management .....	669
Image Options .....	673
Image Backgating Color .....	673
Mask Options .....	674
Image Processing Models .....	674
Plate Options .....	675
Administrator .....	676
Overview .....	676
Data Management Options .....	677
Other Settings .....	679
SAE Configuration .....	680
Rules for enabling the SAE mode .....	682
Enable the SAE mode .....	682

Disable the SAE mode – SAE user .....	684
Disable the SAE mode – local Attune™ administrator .....	686
User Management .....	687
Overview .....	687
Users tab .....	688
System Security tab .....	692
Configuration .....	695
Overview .....	695
Hardware/Virtual laser configuration .....	696
Enable bubble sensor .....	697
Automation communication settings .....	697
Connect device registration .....	698
Unregister a device from the Connect .....	698
Remove individual users from a device on the Connect .....	699
Resources .....	701
Overview .....	701
Access resources .....	701
About Attune™ cytometric software .....	702
Trademarks .....	703
Contact us .....	703
<b>■ CHAPTER 26 User Management .....</b>	<b>704</b>
Access the user management tab .....	705
Add or edit an account .....	705
Add an account .....	705
Edit an account .....	705
Create a user profile .....	706
Edit a user profile .....	708
Set password reminders .....	710
Reset a password .....	711
Change a password .....	712

■	<b>CHAPTER 27</b>	<b>File dialogs</b>	<b>714</b>
	Overview		714
	File dialogs		714
	File Save (Export) dialog		715
	Overview		715
	Dialog behavior		716
	File filters and default names		718
	File Open (Import) dialog		721
	Overview		721
	Dialog behavior		721
	File filters		722
	Folder Browser dialog		723
	Overview		723
	Dialog behavior		724
	File filters and default names		726
■	<b>CHAPTER 28</b>	<b>Dialogs</b>	<b>727</b>
	Overview		727
	Page Setup dialog		728
	Open the dialog		728
	Layout tab		728
	Print dialog		730
	Open the dialog		730
	Print range		730
	Print what options		731
	Print Preview dialog		732
	Open the dialog		732
	Print Preview dialog		733
	Batch Print dialog		734
	Open the dialog		734
	Print dialog options		734
	Print order		735
	Low Disk Space dialog		736
	Export Statistics dialog		737
	Export statistics of selected samples using		737
	Exports statistics of selected views		737
	To		738
	Derived gate dialog		739
	Launch the dialog		739
	Boolean operators		740
	Derived gate dialog		740

Edit gates dialog .....	742
Launch the dialog .....	742
Edit gates dialog .....	743
Confirm file save dialog .....	746
Save As Template dialog .....	747
Open the Save As Template dialog .....	747
Name and Description fields .....	748
Template options .....	748
Template thumbnail .....	749
Template attributes .....	749
Manage templates dialog .....	749
Type filter .....	750
Keyword search .....	750
Template list .....	751
Description .....	752
Thumbnails .....	752
Export .....	754
Import .....	754
Delete .....	755
Plate Mapping dialog .....	756
Plate ID (Barcode) to Template table .....	757
Select All .....	757
Delete and Add buttons .....	758
Import Plates .....	758
Apply Template to Selected Plates .....	758
OK and Cancel buttons .....	759
Select template dialog .....	760
Template keyword search .....	760
Template list .....	761
Template description .....	761
Thumbnails .....	762
Import .....	762
OK and cancel buttons .....	763
Map Sample List Data dialog .....	764
Update sample information dialog .....	766
Create/Edit keyword dialog .....	767
Name and Description .....	768
Type .....	768
Default Value and Decimal places .....	768
Keyword Options .....	769
Experiment Keywords dialog .....	770
Add/Edit .....	770
Remove .....	771
Export Logs for Service dialog .....	772

Select Plates dialog .....	774
Create/Edit plate .....	776
Test Plate dialog .....	778
Run the plate test .....	778
Foil cover present dialog .....	779
Process (Reprocess) Images dialog .....	780
Remove Image Processing Data dialog .....	781
Link Account dialog .....	782
Link account with the QR code .....	783
Link account with the linking code .....	785
Cloud Sign In dialog .....	787
Switch Account dialog .....	790
Connect storage dialog .....	791
Maintenance Required dialog .....	792
System Log dialog .....	793
General tab .....	795
User Log tab .....	798
Maintenance tab .....	800
SAE Signing dialogs .....	804
Overview .....	804
e-Signature workflow .....	804
Request Signature dialog .....	806
e-Signature Request dialog .....	807
Sign Records dialog .....	808
Sign Experiment dialog .....	810
e-Signature Record Report History dialog .....	812
<b>■ CHAPTER 29 Attune™ Database Utility .....</b>	<b>813</b>
Overview .....	813
Startup .....	813
Main application startup .....	813
Startup errors .....	814
Login screen .....	814
Overview .....	814
User sign in .....	815
Login warnings .....	815
Main application .....	816
Overview .....	816
Status tab .....	817
Backup tab .....	818
Restore tab .....	820
Schedule Backup tab .....	822
Reinstall Database .....	823

Database Configuration tab .....	824
Attune™ Database Service .....	825
■ <b>APPENDIX A FCS file reference .....</b>	<b>827</b>
Overview .....	827
FCS HEADER format .....	827
FCS TEXT segment .....	828
Required keywords .....	828
Optional keywords .....	829
Custom keywords .....	831
Storage of user-defined keywords in the FCS file .....	833
Storage of compensation values in the FCS file .....	833
■ <b>APPENDIX B Attune™ Cytometric Software image processing parameters .....</b>	<b>834</b>
Image processing parameters .....	834
Kurtosis .....	836
Skewness .....	836
■ <b>APPENDIX C Data management in Attune™ Cytometric Software .....</b>	<b>837</b>
Overview .....	837
Contents .....	837
Types of data in Attune™ Cytometric Software .....	838
Disk usage .....	839
Data storage warnings .....	841
Change low disk space warning options .....	842
Options for saving user data .....	843
Choosing one or more data archiving methods .....	843
Database backup .....	844
Back up data .....	846
Save and export Plate and Tube Experiment files .....	846
Save and export FCS files .....	848
Save and export FCS and image data .....	850
Save and export images .....	852
Back up data using the Attune™ Database Utility .....	854
Restore data .....	856
Restore data using the Attune™ Database Utility .....	856
Copy a database to a different computer .....	858
Workflow .....	858
Save data using the Attune™ Database Utility .....	858
Copy data and backup file to network or external hard drive .....	860
Install Attune™ Cytometric Software on new computer .....	860
Restore database using the Attune™ Database Utility .....	861

Permanently remove saved data from the Attune™ computer .....	863
Data deletion options .....	863
Delete Tube and Plate Experiment files .....	863
Reinstall a new database to erase existing data .....	866
Delete a user account to remove data .....	868
Permanently remove old baseline and performance test data .....	871
About baseline and performance test data .....	871
Remove baseline and performance test data in Attune™ Cytometric Software version 2.7 or earlier .....	874
Remove baseline and performance test data in Attune™ Cytometric Software version 3.1 and later .....	876
<b>■ APPENDIX D SAE Administrator Console .....</b>	<b>879</b>
SAE module .....	879
Overview of the SAE module .....	879
SAE Administrator Console .....	880
Overview of the SAE Administrator Console .....	880
Open the SAE Administrator Console .....	881
SAE Administrator Console tabs .....	881
Users tab .....	882
Create an SAE user account .....	883
Edit an SAE user account .....	884
Activate a suspended SAE user account .....	884
Disable (inactivate) an SAE user account .....	884
Reset an SAE user password .....	884
Print or view a user report .....	885
Roles tab .....	886
SAE account permissions .....	887
Create a role .....	894
Edit a role .....	895
Delete a role .....	895
Print or view a role report .....	896
System tab .....	897
User Name Settings .....	897
Password Policy .....	898
Account Lockout Policy .....	899
Other Settings .....	900
Audit tab .....	903
Enable Audits .....	904
Audit Settings panel .....	905
Audit Reason Settings panel .....	906
Enable or disable the audit function .....	907
Select items to audit and set the audit mode .....	907
Configure audit reason settings .....	908



e-Signature tab .....	909
Enable or disable the e-Signature function .....	910
Select the actions that require e-Signature .....	910
Configure the meanings of e-Signatures .....	911
Audit History tab .....	913
Action Records audit log .....	914
System Configuration audit log .....	917
Application Objects Records audit log .....	919
Instrument Run Records audit log .....	924
View audit logs (Audit History) .....	924
Settings tab .....	927
Set up SAE messaging notifications .....	928
Configure the SMTP server for email notifications .....	929
Auto archive audit records .....	929
Archive audit records manually .....	930
View or export archived audit records .....	930
Restore archived audit records .....	931
Export configuration .....	932
Import configuration .....	932
Set up the SAE administrator console with application profiles .....	933
Configure user repositories .....	933
User repository settings .....	934
Sign-in with LDAP or federated user repositories .....	934
User repository overview .....	936
■ <b>APPENDIX E Technical reference .....</b>	<b>937</b>
Overton's cumulative statistics .....	937
Kolmogorov-Smirnov test .....	937
Hyperlog™ .....	938
Linear and log color binning reference .....	938
■ <b>APPENDIX F Safety .....</b>	<b>939</b>
Chemical safety .....	940
Biological hazard safety .....	941
<b>Documentation and support .....</b>	<b>942</b>
Customer and technical support .....	942
Limited product warranty .....	942



# About this guide

This user guide describes how to use the Attune™ Cytometric Software to acquire and analyze data using the Attune™ NxT Flow Cytometer or the Attune™ CytPix™ Flow Cytometer and assumes that you have a working knowledge of basic Microsoft™ Windows™ operation.

## Conventions

### Text and keyboard conventions

The following table lists the text and keyboard conventions used in the *Attune™ Cytometric Software User Guide*. For safety alert words and symbols used in this document, see “Safety alert words” on page 30.

Convention	Use
<i>Italics</i>	<i>Italic</i> text highlights new or important terms on their first appearance in the user guide. It is also used for emphasis and for user guide or reference titles. For example: <i>Experiment Explorer</i> lists <i>Experiments</i> in a hierarchal view and functions as an interface for creating new Experiments and recording data.
<b>Bold</b>	<b>Bold</b> text indicates user action. For example: Click <b>Run</b> .
▶	Right arrow symbol (▶) indicates a menu choice and separates successive commands you select from a drop-down or shortcut menu. For example: Select <b>Show Events ▶ All Events</b> .
Ctrl+X	When used with key names, a plus sign means to press two keys simultaneously. For example: Click <b>Ctrl+P</b> .



### Clicking

Unless explicitly stated, clicks are left mouse button clicks. If you have transposed the mouse buttons, the primary click is the left click, even though it can be physically swapped.



## User attention symbols

The following attention symbols are used in the *Attune™ Cytometric Software User Guide*. For safety alert words and symbols used in this document, see “Safety alert words” on page 30.

Symbol™	Use
	<b>Note:</b> Describes important features or instructions, and highlights tips that can save time and prevent difficulties.
	<b>IMPORTANT!</b> Provides information that is necessary for proper instrument operation, accurate installation, or safe use of a chemical.

## Other Attune™ user guides

Guide	Pub. No.
<i>Attune™ Cytometric Software User Guide</i>	MAN0026553
<i>Attune™ NxT Flow Cytometer Quick Reference Guide</i>	100024233
<i>Attune™ NxT Flow Cytometer User Guide</i>	MAN0026547
<i>Attune™ NxT Flow Cytometer Maintenance and Troubleshooting Guide</i>	100024234
<i>Attune™ NxT Flow Cytometer Site Preparation Guide</i>	100024428
<i>Attune™ NxT External Fluid Supply User Guide</i>	100038577
<i>Attune™ NxT External Fluid Supply Quick Reference Guide</i>	100037944
<i>Attune™ NxT Auto Sampler User Guide</i>	100032905
<i>Attune™ CytPix™ Flow Cytometer User Guide</i>	MAN0019440
<i>Attune™ CytPix™ Flow Cytometer Site Preparation Guide</i>	MAN0019443
<i>CytKick™ and CytKick™ Max™ Autosampler User Guide</i>	MAN0018351

Additional resources are available on the Flow Cytometry Technical Resources page at [thermofisher.com/flowresources](http://thermofisher.com/flowresources), where you can find protocols, application notes, and tutorials.



## Safety information

---

**Note:** See “**Appendix E: Safety**” for the complete the chemical or instrument safety information.

---

### Safety alert words

Four safety alert words appear in this user guide at points in the document where you need to be aware of relevant hazards. Each alert word—**IMPORTANT, CAUTION, WARNING, DANGER**—implies a specific level of observation or action, as defined below:

---

**IMPORTANT!** Provides information that is necessary for proper instrument operation, accurate installation, or safe use of a chemical.

---



**CAUTION!** Indicates a potentially hazardous situation that, if not avoided, could result in minor or moderate injury. It is also used to alert against unsafe practices.



**WARNING!** Indicates a potentially hazardous situation that, if not avoided, could result in death or serious injury.



**DANGER!** Indicates an imminently hazardous situation that, if not avoided, will result in death or serious injury. This signal word is to be limited to the most extreme situations.

Except for **IMPORTANT!** safety alerts, each safety alert word in this document appears with an open triangle figure that contains a hazard symbol. These hazard symbols are identical to the hazard symbols that are affixed to the instruments (see “**Symbols on instruments**”).

### SDSs

The Safety Data Sheets (SDSs) for any chemicals supplied by Thermo Fisher Scientific are available to you free 24 hours a day. For instructions on obtaining SDSs, see “**Obtaining SDSs**”.

---

**IMPORTANT!** For the SDSs of chemicals not distributed by Thermo Fisher Scientific contact the chemical manufacturer.

---



# Attune™ Cytometric Software

## Attune™ Cytometric Software overview

The Attune™ Cytometric Software is used to acquire and analyze data using the Attune™ NxT Flow Cytometer or the Attune™ CytPix™ Flow Cytometer and assumes that you have a working knowledge of basic Microsoft™ Windows™ operation.

---

**IMPORTANT!** For instructions on using the Attune™ NxT Flow Cytometer, see the *Attune™ NxT Flow Cytometer User Guide* (Pub. No. MAN0026547) and the *Attune™ NxT Flow Cytometer Quick Reference Guide* (Pub. No. 100024233).

For instructions on using the Attune™ CytPix™ Flow Cytometer, see the *Attune™ CytPix™ Flow Cytometer User Guide* (Pub. No. MAN0019440).

---

The functions of the Attune™ NxT Flow Cytometer and the Attune™ CytPix™ Flow Cytometer are controlled by the Attune™ Cytometric Software. The software is preinstalled to the computer workstation supplied with the instrument, and the Attune™ Software icon (i.e., shortcut) is placed on the computer desktop and under **Start ▶ ThermoFisher ▶ Attune™ Software**.

---

**Note:** To reinstall the Attune™ Cytometric Software for recovery, you must have *Administrator* access (see “Default local Attune™ accounts” on page 42). For instructions to reinstall the Attune™ Cytometric Software, see “Attune™ Cytometric Software installation” on page 33.

---

## About the software

The Attune™ Cytometric Software is a flexible data acquisition and analysis tool that uses a browser view to:

- Design and perform experiments
- Define independent instrument settings and optimize data collection
- Carry out instrument performance checks and track instrument performance
- Acquire and record data
- Manage and process recorded data

## System requirements

A PC workstation running the Attune™ Cytometric Software on a Windows™ 10 64-bit platform is provided with the instrument.

---

**Note:** We recommend that you set the text to small (100%) in the Windows™ display options (96 dpi) for the best display of the software application.

---

## Supported files

- The Attune™ Cytometric Software records and outputs data in FCS (Flow Cytometry Standard) 3.1.
- The Attune™ Cytometric Software Cytometric Software can export data in FCS 3.0 and 3.1 formats.
- FCS formats 3.0 and 3.1 are supported for data analysis.

## Software layout and ribbon bar menu system

### Software layout

The Attune™ Cytometric Software has three main screens, **Login**, **Main Menu**, and **Attune™ Desktop**.

- **Login** – Used for signing into the software. The Login screen is the first screen presented after the main application startup (“SAE module” on page 37).
- **Main Menu** – Used for running and reviewing performance tracking tests, creating new experiments, and accessing the current and stored experiments. Main Menu is the first screen presented after a successful login. For more information, see “Main Menu” on page 46.
- **Attune™ Desktop** – Used for controlling the Attune™ NxT and Attune™ CytPix™ Flow Cytometers to run samples, generate data, and analyze results. **Attune™ Desktop** is the main application window. For more information, see Chapter 3, “Attune™ Desktop”.

### Ribbon bar menu system

The Attune™ Cytometric Software uses the Microsoft™ Ribbon interface. The ribbon bar consolidates related functionality by organizing control elements in logical groups under contextual tabs. Each ribbon tab relates to a type of activity.

The availability of the ribbon tab in the ribbon bar is dependent on the context of the software, i.e., they are contextual tabs that are displayed only when they are needed by the user. Objects that are not available as a function in a specific context are shaded gray in the tab.

For more information, see “Sign in to Cloud” on page 53.

---

**Note:** Only a limited set of ribbon tabs are available on the **Main Menu** (“Main Menu” on page 46) when running Performance Tests (see “Overview” on page 555).

On the **Login** screen (“SAE module” on page 37), only the **File** tab (“File tab” on page 72) is available.

---

## Context menus

The Attune™ Cytometric Software interface also allows the use of *context menus* (also called contextual, shortcut, and pop-up menu), which appear upon user interaction with a right-click mouse operation. A context menu offers a limited set of choices that are available in the current state, or context, of the software and the available choices are actions related to the selected object.

## Keyboard shortcuts

Several keyboard shortcuts are available for use in the Attune™ Cytometric Software.

Keyboard shortcut	Function
Ctrl + N	New Experiment
Ctrl + C	Copy Workspace elements (plots, gates, stats boxes)
Ctrl + V	Paste Workspace elements (plots, gates, stats boxes)
Ctrl + X	Delete Workspace elements (plots, gates, stats boxes)
Ctrl + A	Select all (objects on Workspace)
Ctrl + P	On demand printing
Ctrl + S	Save as template
Ctrl + space bar	Select entire column on Heat map
Shift + space bar	Select entire row on Heat map
Ctrl + mouse wheel	Zoom in/out of Workspace or Overlay view (or galleries)

## Attune™ Cytometric Software installation

The Attune™ Cytometric Software is preinstalled to the computer workstation supplied with the Attune™ NxT and Attune™ CytPix™ Flow Cytometers.

The Attune™ Cytometric Software shortcut is placed on the computer desktop and under **Start ▶ ThermoFisher ▶ Attune™ Software**.

However, if you need to reinstall the software for recovery purposes or install a newer version of the software, follow the instructions below.

---

**Note:** You have the option to install the Attune™ Cytometric Software with or without the Security, Auditing, and e-Signature™ functions (i.e., the “SAE module”), which allow you to configure the software to meet specific requirements for security, audit, and e-Signature™ (21 CFR Part 11) (see “Appendix D, “SAE Administrator Console”).

A license control mechanism in the form of a DESkey device is required for the operation of the Attune™ Cytometric Software after installation. The type of the DESkey device required depends on the installation option selected (i.e., with or without SAE functions) and the number of licenses purchased (single or multiple users).

---

## Guidelines for installation

- **IMPORTANT!** Ensure that all data is backed up to an external storage device before installation.
- The instrument must be powered on and connected to the computer for the firmware updater to run at the end of the installer.
- **DO NOT** update the firmware if the instrument is in a sleep state. The indicator lights on the front of the instrument will fade in/fade out in multiple colors during the sleep state. Power cycle (turn on and off) the instrument before running the firmware updater.
- **DO NOT** launch the software application until all installation steps have been completed.
- **DO NOT** run any other applications while completing these steps.
- No change to existing login credentials will occur during the software upgrade.
- If installing the Attune™ Cytometric Software with the Security, Auditing, and e-Signature™ option, ensure that the computer on which the SAE server is installed is configured with a static IP address. Consult your network administrator for help with checking the IP address configuration.

## Install the Attune™ Cytometric Software

1. Insert the DESkey USB device appropriate for your installation type (i.e., with or without SAE functions) into the computer.
2. Restart or power on the computer and the cytometer.
3. Sign in to Windows™ as:  
User: INSTR-ADMIN  
Password: INSTR-ADMIN

---

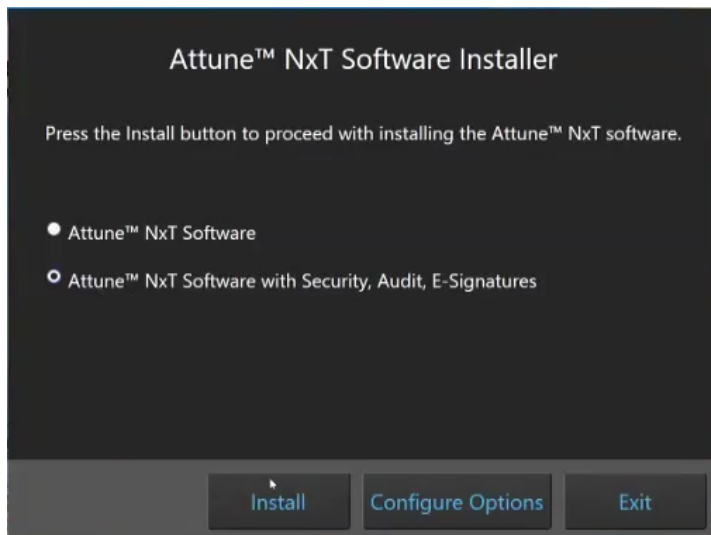
**Note:** This is the default administrator account. If your instrument is on a network, ensure that the administrator privileges have not been removed by your local IT department.

---

4. Unzip (select **extract all files**) the *AttuneNxT\_XX.zip* file to the desktop.  
The XX in the name of the zip file represents the Attune™ Cytometric Software version and must be 5.0 or higher for installation with the Security, Auditing, and e-Signature™ functions.
5. Open the **Attune™ NxT XX** folder, and double-click **SetupAttune.exe** to install the software.



6. In the installer, select **Attune™ Software** (or **Attune™ Software with Security, Audit, E-Signatures**, if SAE is required), then click **Install**.



7. Click **Exit** after the installation is completed successfully.

---

**Note:** When signing into the Attune™ Cytometric Software for the first time after installation, you must sign in as an Administrator. The default username and password for the Administrator are both "admin". After you sign in for the first time, the software will prompt you to change the password.

---

## Update the instrument firmware

After software installation is completed, you must update the instrument firmware. The firmware updater utility launches automatically if the instrument is powered on and connected.

---

**IMPORTANT! DO NOT** update the firmware if the instrument is in a sleep state. The indicator lights on the front of the instrument will fade in/fade out in multiple colors during the sleep state. **Power cycle** (turn on and off) the instrument before running the firmware updater. The firmware update process takes less than 15 minutes.

---

1. On the instrument firmware utility dialog, click **Update Firmware** to update the firmware, then click **OK** to confirm the request to update instrument firmware.
2. After the firmware update is completed, the **Firmware Update Complete** dialog is displayed. Click **Close** to close the dialog to exit the firmware update utility.
3. Cycle the power on the instrument (turn off, then on) to complete the firmware update.
4. Launch the software using login credentials used in earlier software versions.

---

**Note:** If the firmware update is not completed during installation, a prompt to update the firmware is given the first time the application is launched when the instrument is connected and powered up.

---

# 2

## Startup, Login, and Main Menu

### Main application startup

#### Main application startup

To launch the Attune™ Cytometric Software, double-click the **Attune™ Software** shortcut icon on the desktop.

Alternatively, select **Start ▶ ThermoFisher ▶ Attune™ Software**.



---

**IMPORTANT!** A license control mechanism in the form of a DESkey device is required for the operation of the Attune™ Cytometric Software.

---

- If a valid license mechanism is detected when the software is started, the End User License Agreement (EULA) will be displayed on first use of the software (“End user license agreement (EULA)” on page 36).
- If a valid DESkey device is not present, the software displays the appropriate warning message.

#### End user license agreement (EULA)

The End User License Agreement (EULA) is displayed when:

- The Attune™ Cytometric Software is first installed.
- A new version of the Attune™ Cytometric Software is installed.
- A new Windows™ user is attempting to open the Attune™ Cytometric Software.

Upon installation of the Attune™ Cytometric Software, the EULA can also be found with the Licenses directory contained within the application’s installation folder.

Click **I ACCEPT** to accept the terms of the agreement and launch the software.

Click **I DO NOT ACCEPT** to reject the terms of agreement. The EULA dialog is closed and the software is not launched.

## SAE module

The Attune™ Cytometric Software also has an SAE (“Security, Auditing, and Electronic Signature™”) module, which allows you to configure the software to meet specific requirements for security, audit, and e-Signature™ (21 CFR Part 11).

- Use of the SAE mode requires an SAE-specific DESkey device.
- When the SAE mode is enabled (see “SAE Configuration” on page 680), the SAE-specific DESkey is checked upon login when signing into the software as an SAE user.

---

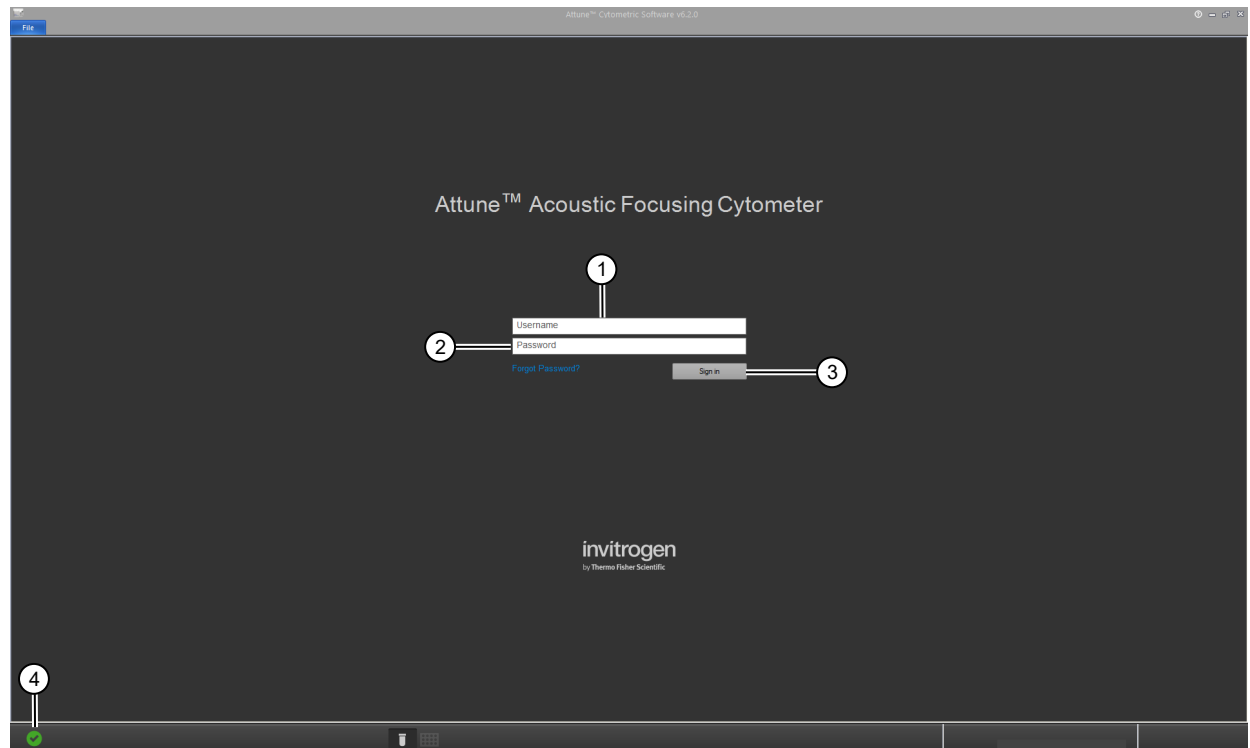
**Note:** For more information on the SAE module and how to configure it to meet specific requirements for security, audit, and e-Signature™, see Appendix D, “SAE Administrator Console”.

---

## Login screen

The **Login** screen is the first screen that is displayed after the **splash** screen when you start the Attune™ Cytometric Software. Each user is required to sign in before being allowed to use the software. The **Login** screen also displays the current Performance Test status.

**Note:** In the SAE (“Security, Auditing, and Electronic Signature”) mode, the sign in button is a split button with a dropdown arrow that enables sign in using the SAE mode or with the local user account (see “Sign in – SAE mode” on page 38).



- ① Username field
- ② Password field
- ③ Sign in button (see “Sign in – standard mode” on page 38)
- ④ Instrument status (see “Instrument status icons” on page 64)

## Sign in – standard mode

1. To sign in the Attune™ Cytometric Software, type a valid username and password in the appropriate text box fields, then click **Sign In**.

Alternatively, you can press the **Enter** key on the computer keyboard after typing the username and password.

2. If the correct username and password are entered, the **Main Menu** is displayed.  
If an invalid username or password is entered, the warning banner displays “Invalid username or password” (see “Invalid username or password” on page 40).

---

**IMPORTANT!** When signing into the Attune™ Cytometric Software for the first time after installation, you need to sign in as an *Administrator*.

The default username and password for the Administrator are both **admin**. For more information about other available account types, see “Default local Attune™ accounts” on page 42.

---

## Sign in – SAE mode

To sign in to the Attune™ Cytometric Software in the SAE (“Security, Auditing, and Electronic Signature”) mode, you need to have an SAE-specific DESkey device installed and the SAE module enabled and configured to allow sign in using SAE credentials. When signing in, you need to have the appropriate SAE user credentials, and meet the SAE account rules for signing in (see “SAE account rules” on page 40).

1. To sign in to the Attune™ Cytometric Software using SAE user credentials, type the valid SAE username and password in the appropriate fields, then click **Sign In**.

Alternatively, you can tap the **Enter** key on the computer keyboard after typing the username and password.

2. To sign in to the SAE-enabled Attune™ Cytometric Software with a local account, click the **drop-down arrow** on the **Sign In** button, then select **Sign in with local account**.

**Note:** If the SAE specific DESkey is not present, you can not sign in and the software displays the “Access Denied” warning dialog.

A screenshot of a login interface. It features two input fields: 'Username' and 'Password'. Below these fields is a blue 'Sign in' button with a small downward-pointing arrow on its right side. At the bottom of the interface is a white button with the text 'Sign in with local account'.

- Click **Yes** to retry the sign in attempt. If the SAE DESkey device is detected, you can sign in as an SAE user.
- Click **No** to close the dialog without signing in as an SAE user.
- If **Yes** is clicked, the dialog persists as long as the SAE Deskey device check fails.

3. To sign in to the Attune™ Cytometric Software in the SAE (“Security, Auditing, and To sign in to the SAE-enabled Attune™ Cytometric Software with a local account, click the **drop-down arrow** on the **Sign In** button, then select **Sign in with local account**.

A screenshot of a login interface, identical to the one above. It features two input fields: 'Username' and 'Password'. Below these fields is a blue 'Sign in' button with a small downward-pointing arrow on its right side. At the bottom of the interface is a white button with the text 'Sign in with local account'.

**IMPORTANT!** When signing into the Attune™ Cytometric Software after the SAE mode is enabled, you need to use the SAE account username and password, and not the Attune™ instrument username and password.

In cases where the SAE server is unavailable or you need service repair of the instrument, you can sign in to the local Attune™ Software account using the **Sign in with local account** dropdown. By default, this option is available if you have an Administrator, System Administrator, or Service account.

To enable local account access to all account types, including User and Advanced User accounts, local Administrator or SAE Administrator must select the **Enable Local Account Access** option in the **Administrator** tab of the **Options** dialog (see “SAE Configuration” on page 680).

Users and Advanced Users are not allowed to login to the local account after SAE is enabled.

Note that the data from local User and Advanced User accounts become inaccessible after the SAE mode is enabled. We recommend that you export all local User and Advanced User data before enabling the SAE mode.

After successful sign in, the following steps are performed:

- If the specified SAE account password has expired, the change password screen is displayed.
- If SAE account is configured to show the password expiration reminder X number of days before the password expires, the password expiration reminder is displayed.
- If the correct username and password are entered, the user login is completed and the **Main Menu** is displayed (see “Main Menu” on page 46).
- If an invalid username or password is entered, the warning banner displays “Invalid username or password” (see “Invalid username or password” on page 40).
- A login audit record is generated (see Note below)
- If it is the first time sign in on the specific Attune™ instrument, a local Attune™ instrument account is created after the successful SAE account sign in. This account is flagged as an SAE user account and is used to allow sign in, if the SAE server is offline.

---

**Note:** In the SAE mode, all successful login events (both local and with SAE credentials) are audited as actions on the SAE server as “Sign In Success” along with the username and role. Local Attune™ user logins are noted with a comment “This is a non-SAE Attune™ user account.” All sign in failures (both local and SAE credentials) are audited as actions on the SAE server as “Sign In Failure” along with the username and role. Local Attune™ user logins are noted with a comment “This is a non-SAE Attune™ user account.”

---

## SAE account rules

Any user with an SAE account can sign in to the instrument if the following conditions are met:

1. User provides valid user credentials.
2. The SAE account is not disabled or suspended by the SAE Administrator.
3. The SAE account is not locked out (due to too many failed sign in attempts).
4. The SAE server is online unless offline sign in has been configured.

## Invalid username or password

- If you enter an invalid username or password, the warning banner on the Login screen displays “Invalid username or password”. You need to then enter the username or the password correctly or the warning banner is displayed again.

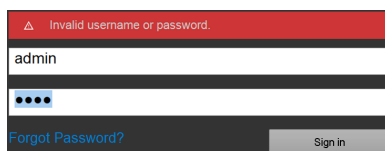


Figure 1 Warning banner for sign in attempt with invalid username or password (local Attune™ account)

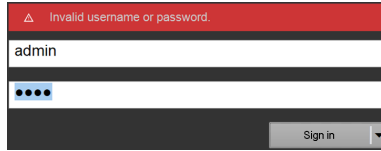


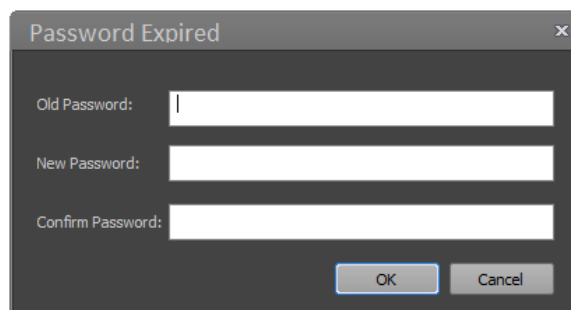
Figure 2 Warning banner for sign in attempt with invalid username or password (SAE account)

- If you attempt to login to an account that has been disabled (i.e., set to inactive) or manually suspended, the warning banner displays “Your account has been disabled. Contact administrator.”

**Note:** SAE user accounts cannot be modified using the Attune™ Cytometric Software (i.e., disabled, deleted, password reset or changed, etc.). All SAE user accounts are managed by the SAE administrator using the **SAE Administrator Console** (see “Roles tab” on page 886” in “Appendix D, “SAE Administrator Console”).

## Expired password

If a password has expired, the Attune™ Cytometric Software prompts you to change your password. You need to enter your old password, your new password, and then confirm the new password.



**Note:** The **Password Expired** dialog is displayed only if you successfully sign in with the correct credentials. If any updates to the password are required, you need to complete them before the Main Menu is displayed.

- If the new password and confirmation password do not match, you are prompted to enter the new password again.
- If the new password that does not meet password complexity requirements, you need to choose another password. The new password must contain at least one uppercase, one lowercase, one numeric, or one non-alpha-numeric character, and must be at least 4 characters long.
- If you enter a password that has been previously reused within the disallowed time frame as specified by the System Administrator, you are prompted to choose another password.
- For more information about user passwords for local Attune™ instrument users, see Chapter 26, “User Management”.

For more information about user passwords for SAE users, see “System tab” on page 897 in Appendix D, “SAE Administrator Console”.

# Account types

## Default local Attune™ accounts

Upon installation, the Attune™ Cytometric Software has three types of default accounts for local Attune™ users, *System Administrator*, *Administrator*, and *Service*.

### System administrator

System Administrator is the highest level account and can define the system security policy for username length, password length, password expiration and lock-out, and the auto lock out time due to system inactivity. The default system administrator account cannot be deleted.

### Administrator

Administrators have full access to the software and can perform a variety of tasks, including managing user accounts, running system diagnostic tests, system decontamination, performance tests and baseline calculations. However, they cannot set system security policy. The default administrator account cannot be deleted.

### Service

Service is a special type of account that gives access to service-only features of the software. The Service™ account cannot be deleted and is only accessible by Thermo Fisher Scientific service team.

---

**Note:** For the permissions assigned to each account type, see “Account permissions” on page 44. For information on how to create additional local Attune™ accounts or delete existing ones, see Chapter 26, “User Management”.

---

## Local Attune™ user accounts

Users can access and modify their own workspace and experiments, and run performance tests, but they cannot run any advanced functions available to other account types. User accounts must be created by the System Administrator or Administrator upon successful installation of the Attune™ Cytometric Software and the initial sign in.

For more information on how to create additional User accounts or delete existing ones, see Chapter 26, “User Management”.



## Default usernames and passwords (local Attune™ accounts)

For the default usernames and passwords for local Attune™ users, see the following table. Note that each account will be forced to reset their password upon first login.

	Default username	Default password
System Administrator	sysadmin	sysadmin
Administrator	admin	admin
User	Set by System Administrator or Administrator when the account is created. <sup>[1]</sup>	Set by default to the username created by System Administrator or Administrator.

<sup>[1]</sup> \*See Chapter 26, “User Management”.

## SAE user accounts

In the **SAE module**, the Attune™ Cytometric Software has four default SAE user account types, **Administrator**, **Advanced User**, **User**, and **Reviewer**. In addition, there is a **No Privileges** role, which is for internal use only by the SAE Administrator Console when setting up user repositories. For more information, see “Roles tab” on page 886).

For the permissions assigned to each default SAE account type, see “SAE account permissions” on page 887.

## Account permissions

For the permissions assigned to each account type, see the following table.

Permission	Description	System Admin.	Admin.	Adv. User	User	SAE User
Advanced Instrument Settings	Enables the adjustment of width threshold, window extensions, and area scaling factor in an Experiment		✓	✓		[1]
Allow Local Login in SAE Mode	Allows a local Attune™ user to login when the SAE mode is enabled	✓	✓	✓	✓	NA
Change Password	Enables a user to change their password	✓	✓	✓	✓	✓
Configure Device	Enables a user to setup a device to connect with Connect Cloud-based platform	✓	✓			[1]
Configure SAE	Allows a user to enable/disable SAE	✓	✓			✓
Create or Edit Global Keyword Access	Enables a user to set and edit global keywords		✓			[1]
Create Plate Definition Access	Enables a user to create and/or duplicate custom plate definitions		✓	✓		[1]
Database Utility Full Access	Enables a user to access the database utility to schedule database and data backups	✓	✓			[1]
Delete Report Column Access	Enables a user to delete plate definitions	✓	✓			[1]
Edit Filters	Enables a user to edit filter configuration filters		✓	✓		[1]
Edit Plate Definition Access	Enables a user to edit plate definitions		✓	✓		[1]
Edit Security Policy	Enables a user to set system security policy (i.e., username and password policies)	✓				[1]

(continued)

Permission	Description	System Admin.	Admin.	Adv. User	User	SAE User
Edit User	Enables a user to modify a user's profile	✓	✓	✓	✓	NA
Export Service Logs	Enables a user to export application logs	✓	✓	✓	✓	✓
Manage User Accounts	Enables a user to create and edit user accounts	✓	✓			[1]
Modify Instrument Configuration	Enables a user to modify instrument configuration		✓	✓		[1]
Run Baseline Calculations	Enables a user to run and reset a baseline		✓	✓		[1]
Run Performance Test	Enables a user to run a performance test		✓	✓	✓	[1]
Run System Decontamination	Enables a user to perform a system decontamination		✓	✓		[1]
Run System Tests	Enables a user to run system diagnostic tests	✓	✓	✓		[1]

[1] SAE user privileges are managed by the SAE administrator using the SAE Administrator Console (see “Roles tab” on page 886” in “Appendix D, “SAE Administrator Console”).

## Main Menu

After you sign in successfully to the Attune™ Cytometric Software, the **Main Menu** is displayed. The **Main Menu** contains the buttons for the most often used commands (**Performance Test**, **New Experiment**, **Import Experiment**, and **Templates**), the **Experiment Explorer** panel, the **Performance test status bar**, and the **Log Out** hyperlink control. If the software is in the Automation mode (“Automation group” on page 94), all buttons and the hyperlink are inactive.



- ① **Main Menu Ribbon Bar** contains the **File** (“File tab” on page 72), **Home** (“Home tab” on page 74), and **Instrument** ribbon tabs (“Instrument tab” on page 88).
- ② **Performance Test** activates the **Performance Test setup** screen in the Performance Test module, which lets you perform the baseline and performance test and view the baseline and performance test reports (see “Overview” on page 555).
- ③ **New Experiment** launches the **New Experiment dialog**, which enables you to create a Plate or Tube-based Experiment, or an Experiment using imported files (see “New Experiment dialog” on page 606).
- ④ **Import Experiment** launches the **File Open (Import) dialog** (“File Open (Import) dialog” on page 721), which enables you to import pre-existing Plate or Tube-based Experiments.
- ⑤ **Templates** launches the **New Experiment from Template dialog**, which enables you to create a new Experiment from a template stored in the database in the account (see “New experiment from template dialog” on page 613).
- ⑥ **Log Out** hyperlink logs out the current user and the software returns to the **Login** screen.
- ⑦ **Instrument status** (see “Instrument status icons” on page 64).

## Log Out

Clicking **Log Out** logs out the current user and the software returns to the **Login** screen (“SAE module” on page 37).

---

**Note:** The **Log Out** link is only enabled when the instrument is not acquiring (see “Acquisition status indicators” on page 344).

---

## Main Menu ribbon bar

The **Main Menu** provides a limited set of options for the user to perform, and contains only the **File** (“File tab” on page 72), **Home** (“Home tab” on page 74), and **Instrument** ribbon tabs (“Instrument tab” on page 88).

- The **Main Menu** ribbon bar is minimized by default. The minimized or expanded state of the **Main Menu** ribbon bar persists on a per user basis.
- When the **Main Menu** ribbon is expanded, the **Home** tab is selected by default.

## Experiment Explorer

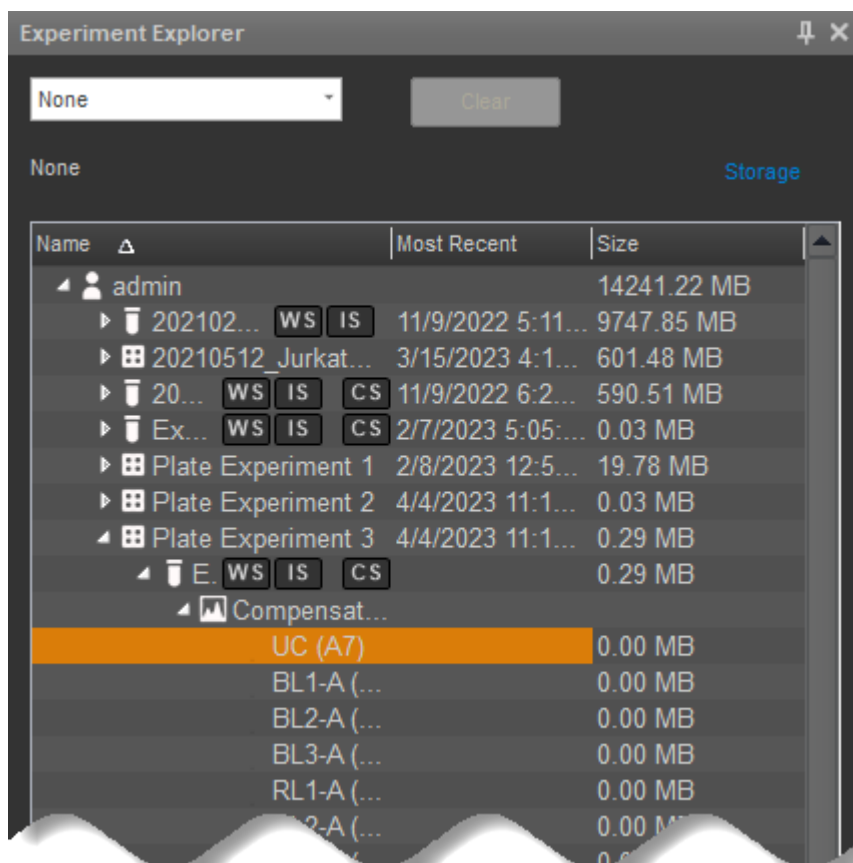
**Experiment Explorer** is used to create, view, and manage experiments, and it is described in detail in Chapter 11, “Experiment Explorer”.

The **Experiment Explorer** panel cannot be undocked or resized in the **Main Menu**.

---

**Note:** **Experiment Explorer** can be undocked and resized in the **Attune™ Desktop**, but not in the **Main Menu**.

---

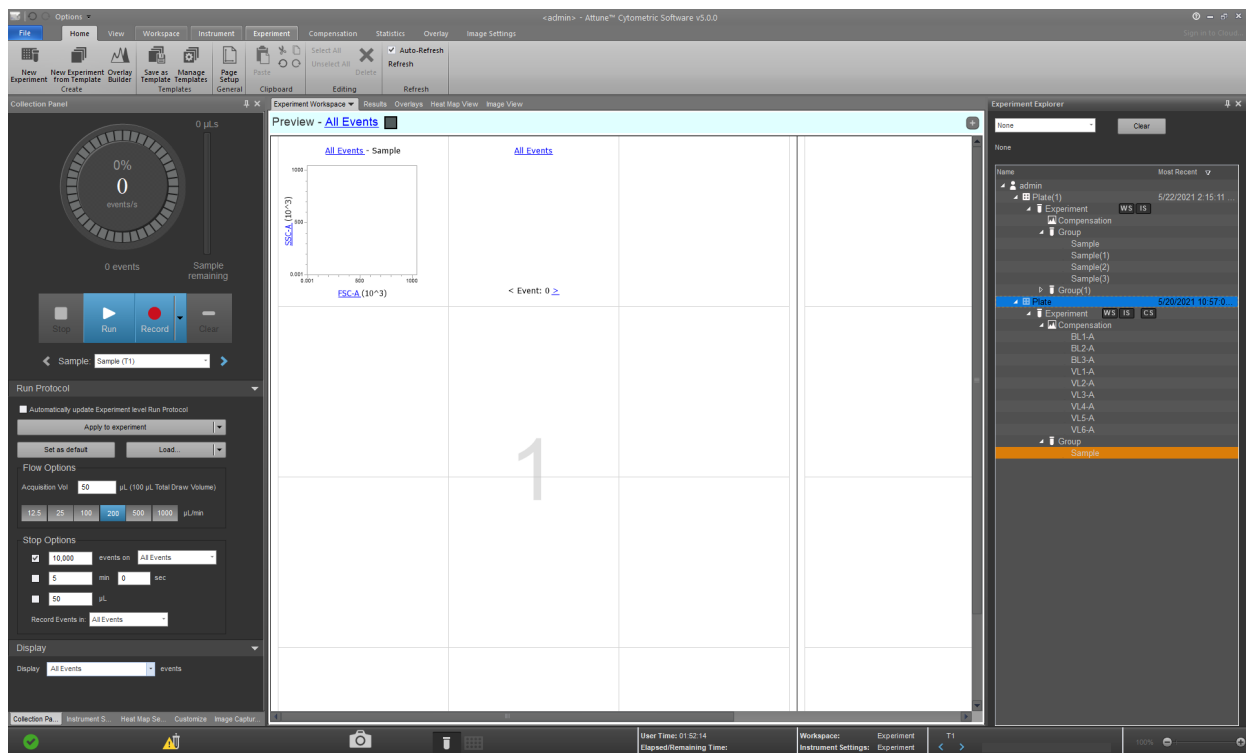




## Attune™ Desktop overview

**Attune™ Desktop** is the main application window, and it is used for controlling the Attune™ NxT and Attune™ CytPix™ Flow Cytometers to run samples, generate data, and analyze results. It consists of the sections listed below.

- **Application title bar** (“Application title bar” on page 51) contains the **Application** button, **Quick Access** toolbar, **Application name**, and **Window Size** buttons.
- **Ribbon bar** (“Sign in to Cloud” on page 53) is organized into a series of tabs that enable access to the main functions of the Attune™ Cytometric Software.
- **Main application workspace** (“Main application workspace” on page 56) accommodates the **Application area**, the adjustable docking panels that contain controls grouped by functionality, the **Experiment Explorer**, and the **Status** notification display.
- **Application status bar** (“Application status bar” on page 63) displays various icons depicting the status of the instrument and the experiment, any alerts relating to the instrument, as well as the progress bar for instrument functions, loading and refreshing. It also contains the **Size slider** (“Size slider” on page 70).



## Attune™ Desktop layout

The layout of the **Attune™ Desktop** uses the ribbon menu system and adjustable docking panels that contain controls grouped by functionality.

The schematic below represents the default layout of the **Attune™ Desktop**, showing the default locations of the various panels and toolbars. Note that the contents of the **Attune™ Desktop** changes depending on the context of the software.

Quick Access Toolbar		Help button	
Labels for Ribbon Bar		Cloud sign in	
Ribbon Bar			
Left panel label	Notification/Status Bar	Right panel label	
Left panel: 1. Collection Panel 2. Instrument Settings 3. Customize Panel 4. Image Capture Settings	Preview	Right panel: Experiment Explorer	
	Tabs for Workspace(s), Results, Heat Map View, Overlays, Sample List, Image View		
	Workspace/Heat Map/Results View		
	Legend (if Heat Map is showing)		
Left panel tabs	Application Status Bar		

---

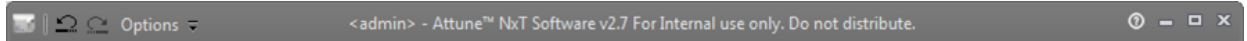
**Note: Application panels** are dockable panels that contain controls grouped by functionality. For more information, see “Application panels” on page 57.


---

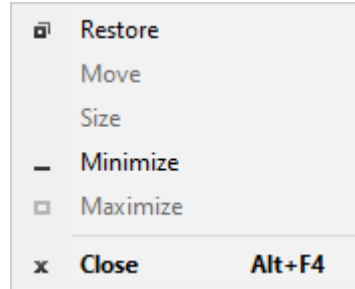


## Application title bar

The **Application title bar** is displayed at the top of the application window, and contains the **Application** button, **Quick Access** toolbar, **Application name**, and **Window Size** buttons.




- **Application** button  opens the standard Microsoft™ Windows™ application size menu.

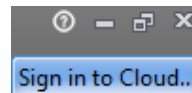


You can also open the application size menu by right-clicking anywhere on the application title bar.

- The **Quick Access** toolbar (“Quick Access toolbar” on page 52) provides easy access to frequently used commands and can be customized to include any commands in the ribbon menu system.



- **Application name** shows the version number of the Attune™ Cytometric Software and the currently logged in username. If these names are too long, they are truncated and ellipses are used.
- **Window size** buttons are the standard buttons used in Microsoft™ Windows™ applications for minimizing, maximizing, and restoring the open application window.
- **Help** () button opens the *Attune™ Cytometric Software User Guide* (Pub. No. [MAN0026553](#)).
- **Cloud sign in** (“Sign in to Cloud” on page 53) button at the lower right of the Application title bar lets you to add the device (instrument or software) to your Connect account, which enables the export and import of files from the Connect cloud-based platform.



## Quick Access toolbar

The **Quick Access** toolbar is a customizable toolbar containing frequently used commands that are independent of the tab currently displayed.








By default, the **Quick Access** toolbar is located above the ribbon bar. However, you can also display the toolbar below the ribbon bar using the **Customize Quick Access Toolbar** dropdown list.

**Note:** Options to customize the **Quick Access** toolbar, to show the toolbar above or below the ribbon, and to minimize the ribbon are also available through the **Ribbon context menu**. For more information, see “Ribbon context menu” on page 55.

### Default buttons

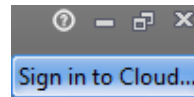
By default, the **Quick Access Toolbar** contains the **Undo** and **Redo** commands, a link to the **Main Menu**, a link to the **Options menu**, and the **Quick Print** command.

Icon	Function
 Undo	Negates the last undoable action performed and brings the application to its previous state. The <b>Undo</b> button is enabled only if there is an undoable action available.
 Redo	Reverses the Undo or advances the application to a more current state. The <b>Redo</b> button is enabled only if there is a redo action available.
 Main Menu	Navigates to the <b>Main Menu</b> (“Main Menu” on page 46). The <b>Main Menu</b> button is enabled only if a user is logged in, and the instrument is not acquiring.
 Options	Opens the <b>Options</b> dialog (Chapter 25, “Options dialogs”), which enables you to customize the Attune™ Cytometric Software by configuring personal settings and changing the default options. The <b>Options</b> button is enabled only if a user is logged in.
	<p>Opens the <b>Customize Quick Access Toolbar</b> dropdown menu.</p> <ul style="list-style-type: none"> <li>• Select the functions displayed on the <b>Customize Quick Access Toolbar</b></li> <li>• <b>Show Quick Access Toolbar Below the Ribbon/Above the Ribbon:</b> Toggles the <b>Quick Access Toolbar</b> to show below or above the Ribbon bar; the text on the menu item toggles between <b>Show Below the Ribbon</b> and <b>Show Above the Ribbon</b> depending on its location.</li> <li>• <b>Minimize the Ribbon:</b> Toggles the Ribbon bar between minimized and restored. When the Ribbon bar is minimized, the tabs are still active. Changing to another tab displays the Ribbon bar.</li> </ul>

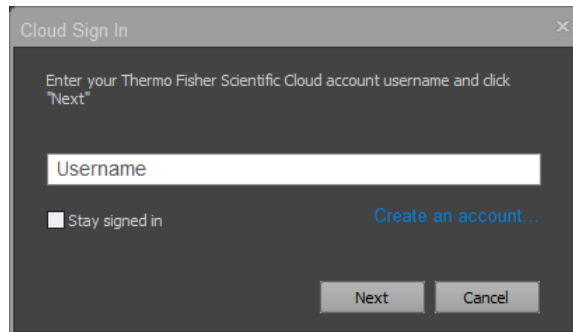
## Sign in to Cloud

**Sign in to Cloud** enables you to connect to the Connect account to export and import of files to and from the Connect cloud-based platform, and to view performance data from a registered device.

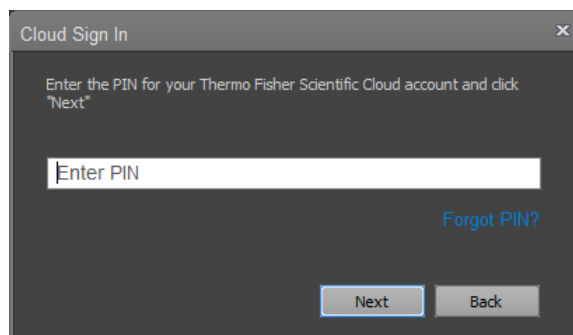
- The **Sign in to Cloud** button appears at the top right of the screen if the device (instrument or software) is registered with the Connect platform. For more information about how to register the instrument or software, see “Connect device registration” on page 698.



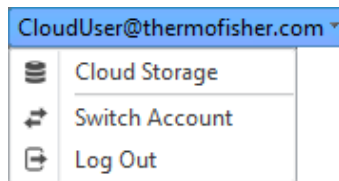
- Click **Sign in to Cloud** to open the **Cloud Sign In** dialog (see “Cloud Sign In dialog” on page 787) and enter the login credentials to sign in to the Connect account.



- If the instrument is registered with the Connect platform but the Attune™ Cytometric Software has not yet been linked with the Connect account, the software displays the **Link Account** dialog (see “Link Account dialog” on page 782).
- If the software and the account are already linked, the **Cloud Sign In** dialog prompts you to enter the four-digit PIN to link the Connect account to the Attune™ Cytometric Software (see “Cloud Sign In dialog” on page 787).



- If you have already signed in to the instrument with a Connect account, the **Sign in to Cloud** button becomes a dropdown button that displays the Connect username on the button text.



- **Cloud Storage:** Opens the **Connect Storage** dialog (see “Connect storage dialog” on page 791), which displays the available storage in the Connect account.
- **Switch Account:** Opens the **Current Attune™ Password Needed** dialog (see “Cloud Sign In dialog” on page 787) to the Connect account password. If the password is correct, the dialog closes and the **Connect Sign in** dialog (“Cloud Sign In dialog” on page 787) opens.
- **Sign out:** Lets you to sign out of the Connect account. When you sign out of the account, the **Remember me** settings in the **Connect Sign in** dialog (“Cloud Sign In dialog” on page 787) are cleared.

## Ribbon bar

The **Ribbon bar** is organized into a series of tabs which represent the main functions of the Attune™ Cytometric Software.

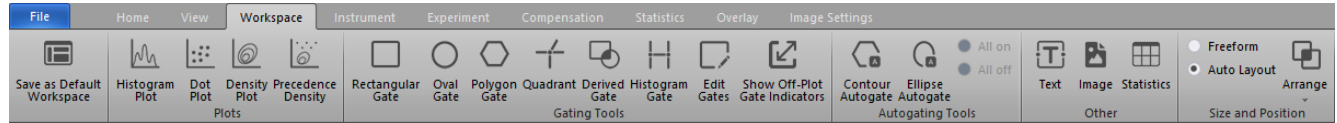


Figure 3 Ribbon bar with the Workspace tab selected (Attune™ NxT Flow Cytometer).

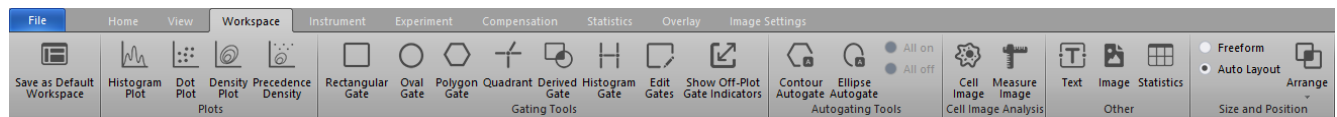


Figure 4 Ribbon bar with the Workspace tab selected (Attune™ CytPix™ Flow Cytometer).

## Ribbon tabs

The availability of the tabs in the ribbon bar is dependent on the context of the software; they are displayed only when the user needs them. Objects that are not available as a function in a specific context are shaded gray in the tab.

Depending on the context of the application and the instrument type (Attune™ NxT or Attune™ CytPix™ Flow Cytometer), the ribbon bar contains one or more of the following tabs:

- **File** tab (page 72)
- **Home** tab (page 74)
- **View** tab (page 77)
- **Workspace** tab (page 81)
- **Instrument** tab (page 88)
- **Experiment** tab (page 95)
- **Compensation** tab (page 99)
- **Statistics** tab (page 103)
- **Overlay** tab (page 110)
- **Image Settings** tab (page 112)
- **SAE** tab (page 114)

**Note:** When performing a Performance Test, only a limited set of tabs are available on the Main Menu. On the **Login** screen, only the **File** tab is available.

**SAE** tab is available only when an SAE user is signed in.

**Image Settings** tab is available only for Attune™ CytPix™ Flow Cytometers.

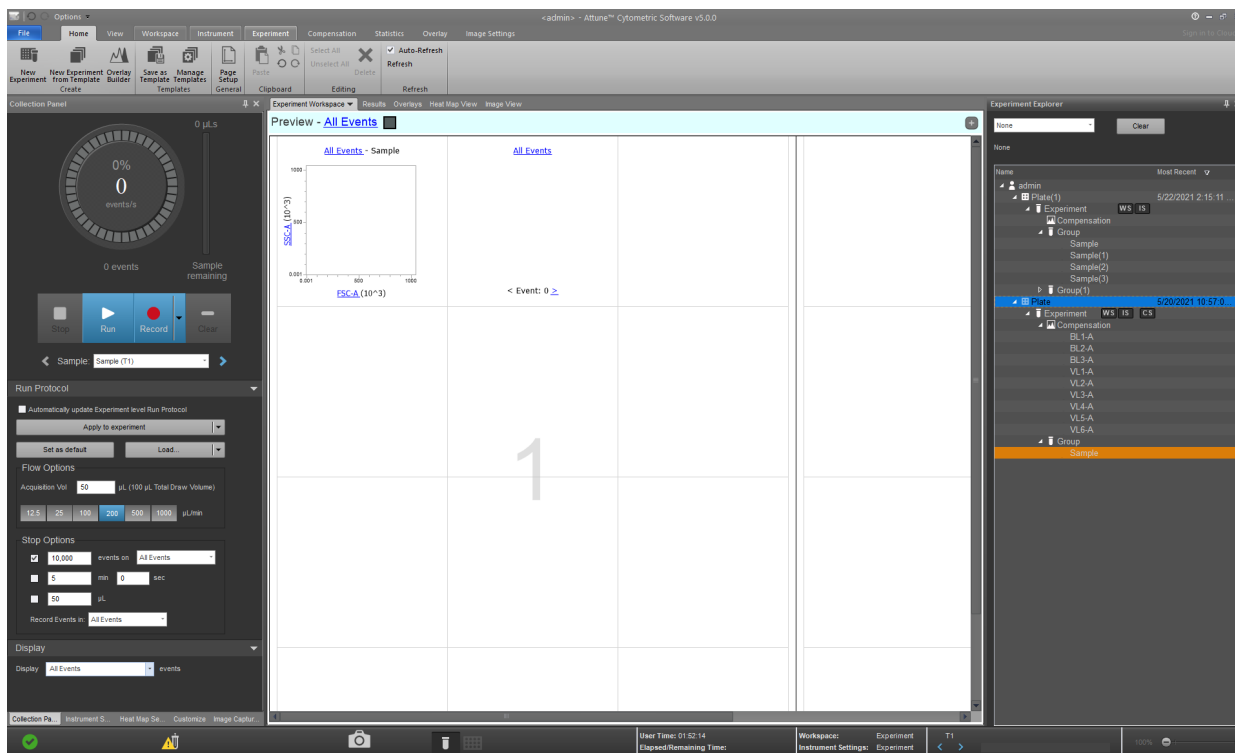
## Ribbon context menu

Right-clicking on a control in the ribbon bar brings up the Ribbon context menu, which contains the options listed below.

- **Show Quick Access Toolbar Below the Ribbon/Above the Ribbon:** Toggles the **Quick Access** toolbar to show below or above the ribbon bar; the text on the menu item toggles between **Show Below the Ribbon** and **Show Above the Ribbon** depending on its location.
- **Minimize the Ribbon:** Checking and un-checking this menu item toggles the ribbon bar between minimized and restored. When the ribbon bar is minimized, the tabs are still active. Changing to another tab displays the ribbon bar.

## Main application workspace

The **Main application workspace** accommodates the **Application area**, the **Application panels**, the **Experiment Explorer**, and the **Status notification** display. The contents of this area change depending on the context of the software.



## Application area

The **Application area** displays the analysis objects (plots, gates, and statistics) that are associated with the current or a saved Sample or Experiment. ext boxes and image files can also be displayed in the **Application area**. The **Application area** contains the following **Application tabs**:

- Workspace view (page 116)
- Heat Map view (page 211)
- Sample List view (page 222)
- Results view (page 228)
- Overlay view (page 234)
- Image view (page 269)

## Application panels

**Application panels** are adjustable docking panels that contain instrument and software controls that are grouped by functionality. The **Main application workspace** contains the following **Application panels**:

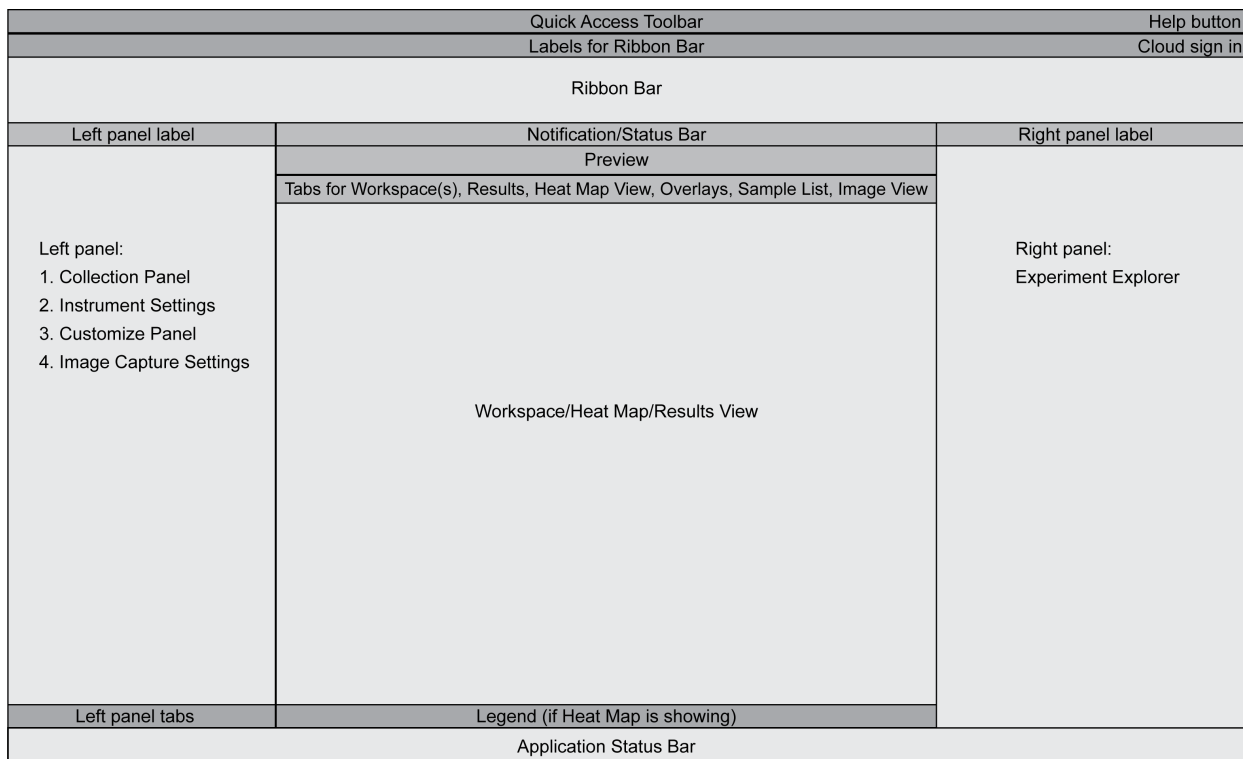
- Experiment Explorer (page 285)
- Collection (page 340)
- Instrument Settings (page 382)
- Customize (page 420)
- Image Capture Settings (page 491)
- FCS Information (page 503)

---

**Note:** The **Application** tabs and the **Application panels** are displayed in the **Main application** workspace as described above, except in the following application-specific modules:

- **Login screen** – All panels are hidden.
  - **Main Menu** – All panels are hidden except for Experiment Explorer, which cannot be undocked or resized.
  - **Performance Test/Baseline Test** – All panels are hidden.
  - **Image View** tab and **Image Capture Settings** panel are available only when using an Attune™ CytPix™ Flow Cytometer or when an Experiment that was created with an Attune™ CytPix™ Flow Cytometer is active.
-

## Docking locations



- By default, the **Collection**, **Instrument Settings**, and **Customize** panels are docked to the left of the **Application area** in the order listed.
- The **Experiment Explorer** is by default docked to the right of the **Application area** (see the schematic above).
- The **FCS Information** panel is set as a floating panel by default (i.e., with its “Dockable” attribute deselected; see “Title bar context menu” on page 62).
- The **Workspace**, **Results**, **Heat Map View**, **Overlays**, **Sample List**, and **Image View** tabs are docked to the **Application area**. By default, they appear in the order listed when docked.

---

**Note:** **Image View** tab and **Image Capture Settings** panel are available only when using an Attune™ CytPix™ Flow Cytometer or when an Experiment that was created with an Attune™ CytPix™ Flow Cytometer is active.

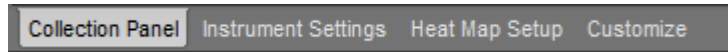
---

- You can reorder all tabs except the **Workspace** tab, which always appears first in the list. To reorder the tabs, click a tab and drag it to different position.
- When multiple panels are docked together to the left or the right of the **Application area**, the name of the currently displayed panel appears in the **container title bar** (i.e., the title bar immediately above the panel). When docked to the **Application area**, no **title bar** is displayed.
- When docked to the **Application area**, tabs for each docked panel appear on the top edge of the **Application area**.

When docked to the left or right of the **Application area**, docked panels are displayed as tabs on the bottom of the panel, unless only one panel is in the **panel container**.



- The tab containing the currently open panel is highlighted.



- **Application panels** that are docked to the **Application area** are shown as tabs.
- The layout of dockable panels persists for each user.

## Docking and undocking panels

1. To undock an **Application panel**, click and drag the title bar of the panel container. While dragging, docking icons are displayed in the left, center, and right of the **Application area** to show where the panel can be docked.

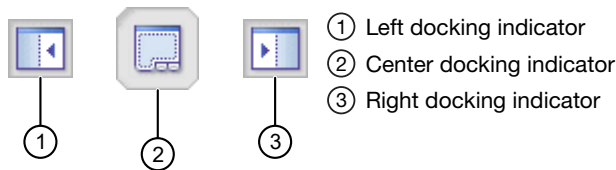


Figure 5 Application panel docking indicators

2. When you place the pointer over one of the **docking indicators**, the docking location of the panel is indicated in light blue (in the example below, to the right of the **Application area**).

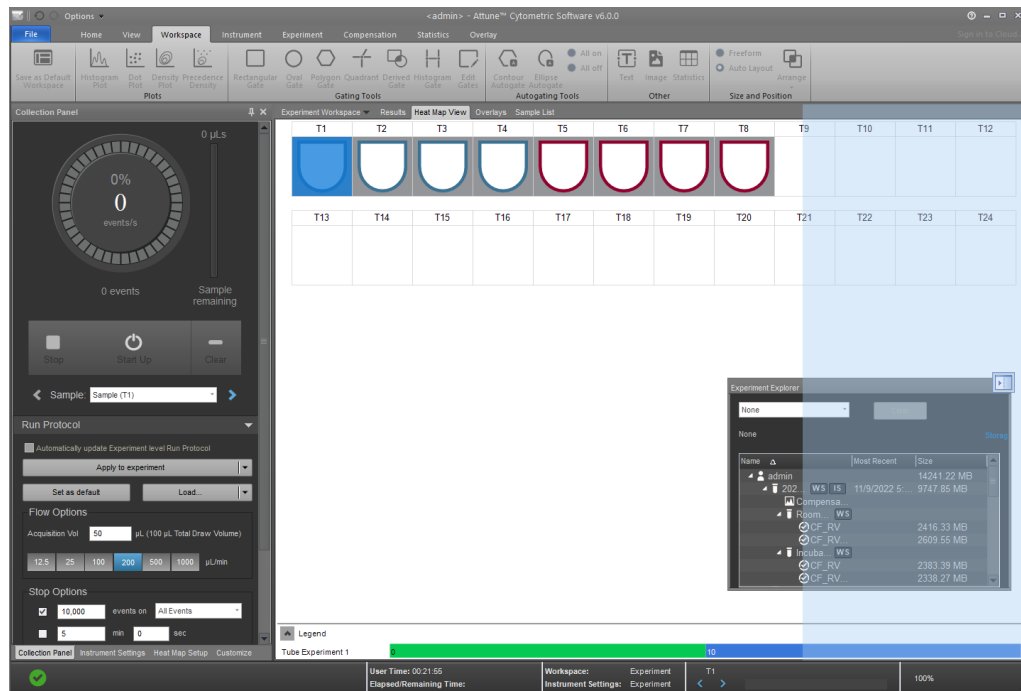


Figure 6 Application panel docking location

3. Release the panel on the desired docking indicator to dock it to the area indicated.

## Docking and undocking tabs

1. To undock an **Application tab**, select and drag the individual tab.
2. Hover over an existing panel to show the **docking indicator**, which enables you to dock the selected tab to any of four sides of the panel (if allowed) or to add it as a tab to the full-size panel.



3. Release the tab on the desired docking icon to dock it to the left or right of the **Main application workspace** or to any of the four docking points of another panel.

You can also drop a selected panel onto the **title bar** of an existing panel to dock it as a full-size panel, which is then added as a tab.

---

**Note:** **Heat Map** and **Results** tabs can be docked anywhere on the **Main application workspace**. All other tabs and panels are limited to left and right only docking rules.

---



## Floating panels

- Double-click the **title bar** of a docked panel or tab to float the panel to its default position.
- A panel that has been docked from a floating position, returns to the same floating position when undocked.
- If multiple floating panels are docked together before docking the entire container to the left or right of the **Main application workspace**, clicking the **title bar** of the docked container refloats the container in its entirety to its previous position.
- Double-click a floating container **title bar** or **panel tab** to redock the container or panel to its last docked position.

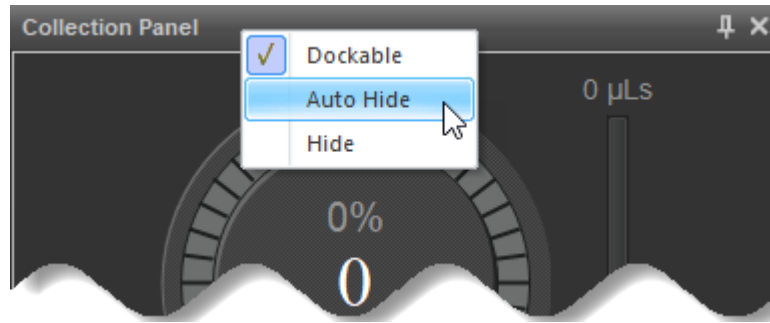
## Auto Hide

**Auto Hide** option reduces a panel to a tab icon when the focus is lost by clicking anywhere outside the menu.

**Push Pin** icon is used for turning the **Auto Hide** option on and off, and it is available only when a panel is docked. The **Push Pin** icon changes to show the current state of the panel.

-  Panel pinned – Auto Hide is turned off.
-  Panel unpinned – Auto Hide is turned on

Alternatively, you can turn the **Auto Hide** option on and off using the **Title bar context menu** (see “Title bar context menu” on page 62).



- When hidden, the panel name appears as a tab on the side of the **Main application workspace** where the panel was docked.
- Clicking on a tab only displays the selected panel.
- Clicking on the **Push Pin** of an open panel restores all panels with which it was docked.

## Close

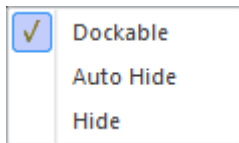
Use the **Close** icon (X) to close the current panel. This deselects the current panel in the **Show Panels** group of the **View ribbon tab** (see “Show Panels group” on page 78).

Alternatively, you can select the **Hide** option using the **Title bar context menu**.

- The only way to redisplay a closed panel is to reselect the relevant checkbox in the **Show Panels** group of the **View ribbon bar** (“Show Panels group” on page 78).
- If multiple panels are docked together and the panel is docked, the **Close** icon only hides the current panel. If the panel is undocked the entire panel is closed.

## Title bar context menu

Right-clicking on the title bar of a panel displays the **Title bar context menu**, which contains the options listed below.



The current state of the panel is indicated in the context menu by a check mark.

- **Dockable:** Enables the docking of a panel. A panel cannot be docked unless this option is checked.

Unchecking the **Dockable** option for a docked panel floats that panel, which cannot be redocked until the **Dockable** option is reselected.

**Dockable** option is not available when the **Auto Hide** option is selected, because a panel must be docked to enable auto-hiding.

- **Auto Hide:** Reduces a panel to a tab icon when the focus is lost by clicking anywhere outside the menu. This has the same action as selecting the **Push Pin** icon in the panel header (see “Auto Hide” on page 60).

The **Auto Hide** option is only available when the panel is docked, not floating.

- **Hide:** Hides the current panel. When hidden, the panel name appears as a tab on the side of the **Main application workspace** where the panel was docked.

Selecting the **Hide** option from the context menu deselects the current panel in the **Show Panels** group on the **View tab** of the ribbon bar (see “Show Panels group” on page 78).

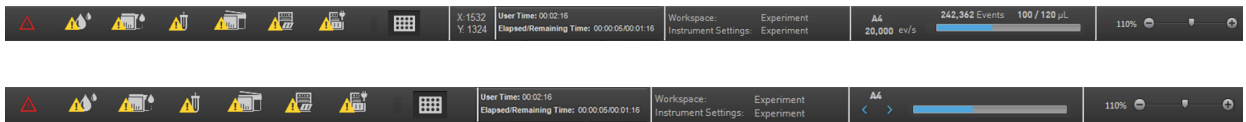
The only way to redisplay a hidden panel is to reselect the relevant checkbox in the **Show Panels** group on the **View tab** of the ribbon bar.

If multiple panels are docked together, the **Hide** option only hides the current panel.

## Application status bar

The **Application status** bar indicates the status of the cytometer and its peripherals and provides feedback on the system using indicator icons. It also contains additional controls to navigate between samples.

The configuration of the **Application status** bar depends on whether the instrument is acquiring (top) or whether it used for analysis (bottom).



The **Application status** bar is described in parts as listed below:

- **Instrument status and alerts** (page 63)
- **Camera status** (page 66)
- **Connect status** (page 66)
- **Tube and Plate modes** (page 66)
- **Time** (page 68)
- **Coordinates** (page 68)
- **Workspace and Instrument Settings** (page 69)
- **Acquisition** (page 69)
- **Analysis** (page 69)
- **Size slider** (page 70)

---

**Note:** Clicking indicator icons does not affect the instrument. They are only used for providing feedback on the status of the system.

---




## Instrument status and alerts

The image below shows the fixed positions of **Instrument Status** and **Alerts** indicator icons on the status bar. If a specific indicator icon is not displayed, then a gap is left in its position.




Note that some indicator icons that are displayed depend on the instrument model and system configuration (for example, camera connected/disconnected icon is available only when an Attune™ CytPix™ instrument is in use).






## Instrument status icons

Icon	Status
 <b>Instrument Idle</b>	<p>The <b>Instrument Idle</b> icon is displayed when the instrument is connected and there are no alarms, or if an instrument has never been connected.</p> <p>The tooltip displays "Idle".</p>
 <b>Instrument Busy</b>	<p>The <b>Instrument Busy</b> icon is displayed when the instrument is connected and it is not idle.</p> <p>The tooltip displays "Instrument busy".</p>
 <b>Alarm</b>	<p>The <b>Alarm</b> icon is displayed when there is an instrument error state.</p> <p>The tooltip displays the relevant error message. If multiple error messages are present, each error message is separated by a comma.</p>



## Instrument alert icons

Icon	Status
 <b>Instrument Fluid</b>	<p>The <b>Instrument Fluid</b> icon is displayed when a fluid level warning is given.</p> <p>The tooltip can contain more than one message concerning the fluidics, using commas between messages:</p> <ul style="list-style-type: none"> <li>• "Shutdown solution tank low"</li> <li>• "Wash tank low"</li> <li>• "Focusing fluid tank low"</li> <li>• "Waste tank full"</li> <li>• "Auto sampler focusing fluid tank low"</li> <li>• "Auto sampler waste tank full"</li> </ul>
 <b>Plate Leak Detected</b>	<p>The <b>Plate Leak Detected</b> icon is displayed when a leak in the Auto Sampler has been detected.</p> <p>Tooltip displays "Auto Sampler leak detected".</p>
 <b>Instrument Leak Detected</b>	<p>The <b>Instrument Leak Detected</b> icon is displayed when a leak in the instrument has been detected. Tooltip displays "Instrument Leak Detected".</p> <p>If a leak is also detected in the Auto Sampler, the Tooltip displays "Instrument Leak and Auto Sampler Leak Detected".</p>

(continued)


Icon	Status
 <b>Tube Lifter Down</b>	<p>The <b>Instrument Tube Lifter</b> icon is displayed when the tube lift is not in the up position.</p> <p>If the tube lifter is down, the Tooltip displays “Tube lifter is down”.</p> <p>If the tube lifter is in the intermediate state, the Tooltip displays “Tube lifter is intermediate”.</p>
 <b>Instrument Cover Open</b>	<p>The <b>Instrument Top Cover</b> icon is displayed when the top cover of the cytometer is not properly closed.</p> <p>Tooltip displays “Top cover is open”.</p>
 <b>Auto Sampler Door Open</b>	<p>The <b>Auto Sampler Door Open</b> icon is displayed when the Auto Sampler door is not properly closed.</p> <p>Tooltip displays “Auto Sampler door is open”.</p>

### Connection status icons

Icon	Status
 <b>Cytometer Not Connected</b>	<p>The <b>Cytometer Not Connected</b> icon is displayed if the cytometer is not connected.</p> <p>The tooltip displays “Attune™ not connected”.</p> <p>If both the cytometer and the Auto Sampler are not connected, the Tooltip displays “Attune™ and Auto Sampler are not connected”.</p>
 <b>Auto Sampler Not Connected</b>	<p>The <b>Auto Sampler Not Connected</b> icon is displayed if the Auto Sampler is not connected.</p> <p>The Tooltip displays “Auto Sampler is not connected”.</p>

**Note:** The connection status icons are only displayed if the software has previously been connected to an instrument as indicated by the database flag, and the conditions described above are met.



### Automation mode icon

Icon	Status
 <b>Automation Mode</b>	<p>The <b>Automation Mode</b> icon is displayed when connected to an instrument and Automation Mode is enabled (“Automation group” on page 94).</p> <p>The Tooltip displays “Automation mode”.</p>

:




## Camera status

The Camera status icons are displayed only if an Attune™ CytPix™ Flow Cytometer is in use.

Icon	Status
 Camera connected	If the Attune™ CytPix™ camera is connected and functioning properly, the <b>Camera connected</b> icon is displayed.
 Camera is not connected	If the Attune™ CytPix™ camera is not functioning properly or the SBC network connection has failed, the <b>Camera is not connected</b> icon is displayed.


## Connect status

The Connect status icons are displayed only if the software or the instrument has been registered to a Connect cloud-based platform account.

Icon	Status
 Registered	If the instrument is registered, the <b>Connect</b> icon is displayed.
 Registered and Signed in	If the instrument is registered and the user is signed in to their Connect account, the <b>Connect</b> icon is displayed with the user icon.
 Connect connection error	If the instrument is registered with a Connect account, but the connection cannot be established, the <b>Connect</b> icon is displayed with the warning icon to indicate a connection error.

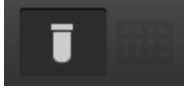


## Tube and Plate modes

When connected to an instrument, either the **Tube** or the **Plate** icon is always displayed.

Icon	Status
 Tube mode – Plate present	When the Experiment is set to Tube mode and the Auto Sampler contains a plate, both the Tube and Plate icons are displayed and the Tube icon is selected. The Tooltip displays “Tube mode”.

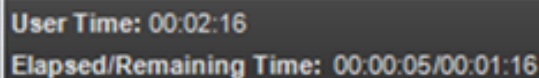


(continued)

Icon	Status
 <p data-bbox="269 426 488 485"><b>Tube mode – Plate absent</b></p>	<p data-bbox="534 321 1485 384">When the Experiment is set to Tube mode and the Auto Sampler does not contain a plate, the Tube icon is displayed and the Plate icon is greyed out.</p> <p data-bbox="534 401 919 432">The Tooltip displays “Tube mode”.</p>
 <p data-bbox="269 621 488 680"><b>Plate mode – Plate present</b></p>	<p data-bbox="534 516 1458 579">When the Experiment is set to Plate mode and the Auto Sampler contains a plate, the Plate icon is selected and the Tube icon is greyed out.</p> <p data-bbox="534 596 1271 627">The Tooltip displays “Plate mode, plate present in Auto Sampler”.</p>
 <p data-bbox="269 816 488 875"><b>Plate mode – Plate absent</b></p>	<p data-bbox="534 711 1485 774">When the Experiment is set to Plate mode and the Auto Sampler does not contain a plate, the Plate icon is selected, but it is greyed out.</p> <p data-bbox="534 791 1256 823">The Tooltip displays “Plate mode, no plate in the Auto Sampler”.</p>

## Time

The **Time** panel displays the **User Time** and **Elapsed/Remaining Time**.



User Time: 00:02:16  
Elapsed/Remaining Time: 00:00:05/00:01:16

- **User Time:** Shows the length a user has been logged on to the instrument in the HH:MM:SS format. The time resets after the user logs out.
- **Elapsed/Remaining Time:** When an Experiment (Plate, Tube, or Manual well) is processing (Run or Record), the status bar displays the **Elapsed/Remaining Time** in the HH:MM:SS format. The elapsed time counter only increments while processing. If the Experiment is paused, the elapsed time counter continues to increase. If the Experiment is paused, the remaining time counter also pauses. When the Experiment stops, the **Elapsed/Remaining Time** also stops and the time persists in the status bar until another run is initiated (Tube or Plate, Run or Record) or the active Experiment is changed (this includes going to another application view such as the **Main Menu**, **Login**, etc.). When the counter is reset or cleared, the time becomes blank.

---

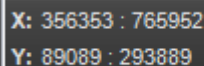
**Note:** If the timer reaches 99:59:59 for the **User Time** or the **Elapsed/Remaining Time**, the displayed timer stops at this value. However, the time recorded in the log file displays the correct time.

---

- The Run Protocol times for the **Remaining Time** counter are calculated as described below:
  - If the stop criteria are based on Event, the estimated **Remaining Time** defaults to the maximum possible Run time for the Normal well, Tube, or Manual well, given the specified Acquisition volume and Flow rate.
  - For stop on Volume, the time is calculated based on the time it takes to process the given volume at the specified Flow rate.
  - For stop on Time, the time for a given Normal well, Tube, or Manual well is based on the specified Stop time.
  - When processing a Normal well, Tube, or Manual well, the estimated time remaining is decremented based on the current rate towards completion of the Well.
  - For Plate experiments, extra time is added to the estimate of each well based on the number of rinses, number of mixes, Wait before record options, Mix mode, Aspiration rate, and Screening mode parameters (these time constants are model dependent).
  - For Plate experiments, the estimated remaining time is recalculated after each well.

## Coordinates

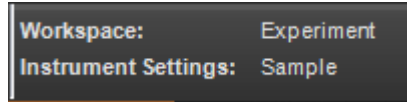
The **Coordinates** show the X and Y coordinates of the selected gate as described in “Move and resize gates” on page 148.



X: 356353 : 765952  
Y: 89089 : 293889

## Workspace and Instrument Settings

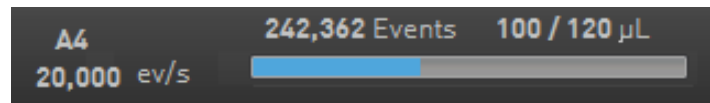
The **Workspace and Instrument Settings** panel displays the open Workspace and the Instrument settings for the active Sample.

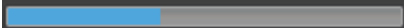


- **Workspace:** Shows the Experiment-level Workspace name during acquisition. Depending on the Workspace selected from the **Workspace** dropdown (“Workspace behavior” on page 117) for the active Sample, this is **Experiment**, **Group**, or **Sample**.
- **Instrument Settings:** Shows the lowest level Instrument Settings, either **Experiment** or **Sample**.

## Acquisition status

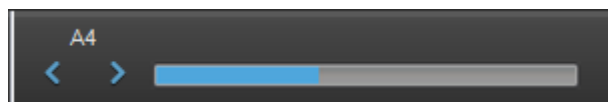
The **Acquisition status** is visible only during acquisition and provides information about the following parameters:



- Well location for a Plate or the Tube ID **A4**
- Calculated event rate (number of events/acquisition time) **20,000 ev/s**
- Number of events acquired at the current time **242,362 Events**
- Current acquired volume and the total available volume **100 / 120 µL**
- Acquisition progress bar showing the progress towards the *Stop condition*, where 100% activates the stop condition 

## Analysis status

The **Analysis status** panel is visible only when an active Experiment is open for analysis and the instrument is not acquiring. It consists of **Sample location (Well location or Tube ID)**, **navigation buttons**, and the **Analysis progress bar**.



- **Well location** for a Plate or the **Tube ID** is displayed for active Experiments.
- The navigation buttons allow navigation between Samples for the current Experiment.
- The **Left navigation** button is disabled when the first Sample in an Experiment is active.
- If the last Tube Sample in an Experiment is active, and the 400 Sample limit has not been reached, the **Right navigation** button is active and adds a new Tube Sample, if clicked. The Tooltip for the button displays “Create new Sample”.

In all other cases when the navigation buttons are enabled, the Tooltip displays either the previous or the next sample name as appropriate.

- If the 400 Sample limit has been reached, the **Right navigation** button is disabled.
- The **Analysis progress bar** is displayed during instrument functions and when loading or refreshing data.
- If multiple operations requiring a progress bar occur simultaneously, the **Analysis progress bar** displays the operation requiring the longest time to be completed.
- In the collapsed state, the priority for display is: instrument function, refreshing, loading. As operations are completed, remaining operations are brought to the foreground.
- The **Analysis Progress bar** is not displayed for tasks that take less than 3 seconds to complete.

## Size slider

The **Size** slider adjusts the size of a selected panel in the **Main Application Window**.



- The **Size** slider is contextual for the current view. If a panel that is not zoomable is in focus, the **Size** slider is disabled. The views for which this is applicable are:
  - **Workspace**
  - **Preview**
- The zoom settings persist per Experiment.
- Size adjustments apply only to the screen display and do not affect the print view.
- The scale of the slider runs from 10% to 400% magnification, with tick marks shown at 10%, 100%, and 400% magnification.
- The scale shows the current scale value rounded to the nearest percent.
- Click the + button to increase the magnification to the nearest whole 10%. For example, from 93% to 100%.
- Click the – button to decrease the magnification to the nearest whole 10%. For example, from 93% to 90%.
- Drag the **Slider** control to adjust the plot size according to its position. The minimum drag rate is a single percent.
- A left-click on the scale, lower to upper tick, sets the **Slider** control to match the horizontal position.

---

**Note:** You can also adjust the **Size** slider by holding the **Ctrl** key down and turning the mouse wheel. This adjusts the view that is currently in focus or underneath the mouse cursor at the time you make the adjustment.

---

# 4

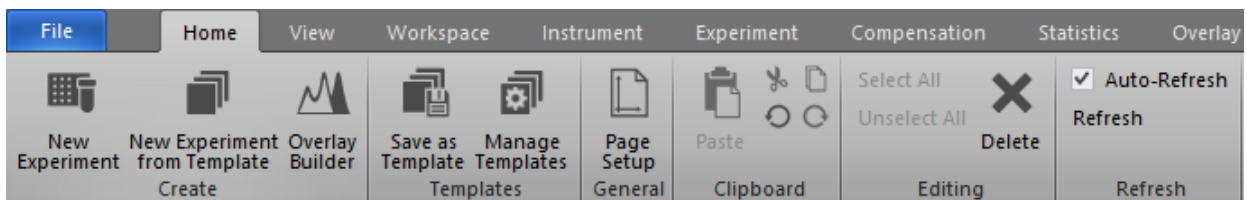
## Ribbon tabs

### Overview

#### Ribbon tabs

**Ribbon tabs** are contextual tabs that combine related software and instrument functionality in the ribbon bar.

The availability of the ribbon tabs in the ribbon bar is dependent on the context of the software, i.e., the tabs are displayed only when they are needed. Similarly, objects and controls that are not available as a function in a specific context are shaded gray in the tab.



Depending on the context of the application and the instrument type (Attune™ NxT Flow Cytometer or Attune™ CytPix™ Flow Cytometer), the ribbon bar contains one or more tabs:

- **File** (page 72)
- **Home** (page 74)
- **View** (page 77)
- **Workspace** (page 81)
- **Instrument** (page 88)
- **Experiment** (page 95)
- **Compensation** (page 99)
- **Statistics** (page 103)
- **Overlay** (page 110)
- **Image Settings** (page 112)
- **SAE** (page 114)

---

**Note:** When performing a Performance Test, only a limited set of tabs are available on the **Main Menu**.

---

On the **Login** screen, only the **File** tab is available.

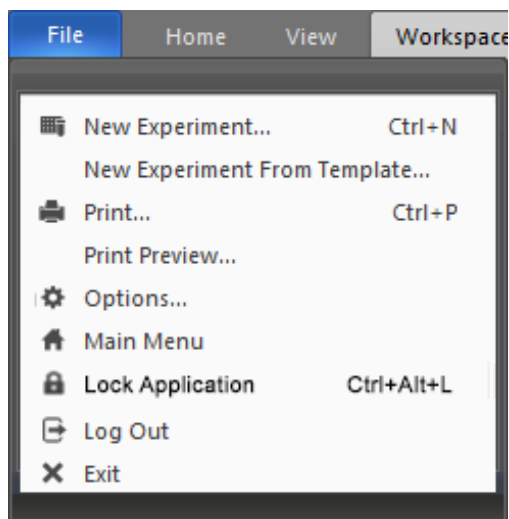
**SAE** tab is available only when an SAE user is signed in.

The **Image Settings** tab is available only for the Attune™ CytPix™ Flow Cytometer or when an Experiment that was created with an Attune™ CytPix™ Flow Cytometer is active.

## File tab

### File menu

**File** tab is the only ribbon bar tab that opens a dropdown menu.



**File** tab opens the **File** dropdown menu, which provides these options:

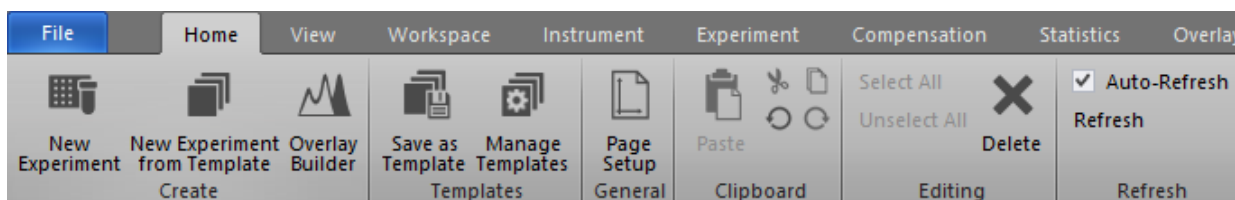
- **New Experiment:** Opens the **New Experiment** dialog (“New Experiment dialog” on page 606), which enables you to create a Tube Experiment, a Plate Experiment, or an Experiment using imported files. The **New Experiment** option is enabled only when the instrument is not acquiring.
- **New Experiment from Template:** Opens the **New Experiment from Template** dialog (“New experiment from template dialog” on page 613), which enables you to create a new Experiment from an existing template with predefined settings. This option is enabled only when the instrument is not acquiring.
- **Print:** Opens the **Print** dialog (“Print dialog” on page 730), which is a modified Windows™ **Print** dialog. It contains the standard Windows™ print commands, as well as the option to select **Workspace, Heat Map, Results, Overlays, and Reports** for printing. The **Print** option is enabled only when the instrument is not acquiring. It is not enabled on the **Main Menu** (“Main Menu” on page 46).
- **Print Preview:** Opens the **Print Preview** dialog (“Print Preview dialog” on page 733), which displays the document as it will appear when printed and contains the controls for various printing options. The **Print Preview** option is enabled only when the instrument is not acquiring.
- **Options:** Opens the **Options** dialog (Chapter 25, “Options dialogs”), which is used to customize the Attune™ Cytometric Software. Some options are user-specific, while others are application-specific (i.e., global to all users) and customizable only by an authorized user. This option is enabled only when the instrument is not acquiring.
- **Main Menu:** Opens the starting **Main Menu** (“Main Menu” on page 46). This option is enabled only when the instrument is not acquiring.
- **Lock Application:** Locks the Attune™ Cytometric Software. This option is visible only if the software is configured to use the **SAE mode**. This option is enabled when a user is signed in. If no user is signed in, the command is disabled.

- **Log Out:** Closes the Attune™ Cytometric Software and logs out the current user, then displays the **Login** screen (“SAE module” on page 37). During Log Out, all FCS files that have not been backed up are backed up automatically. The **Log Out** option is enabled only when the instrument is not acquiring.
- **Exit:** Closes the Attune™ Cytometric Software. This option is enabled only if the instrument is idle.

## Home tab

**Home** tab provides Experiment setup and Output options, and it is organized into six functional groups:

- **Create** (“Create group” on page 74)
- **Templates** (“Templates group” on page 74)
- **General** (“General group” on page 75)
- **Clipboard** (“Clipboard group” on page 75)
- **Editing** (“Editing group” on page 76)
- **Refresh** (“Refresh group” on page 76)



### Create group

**Create** group enables you to create a new Experiment.

**New Experiment:** Opens the **New Experiment** dialog (“New Experiment dialog” on page 606), which enables you to create a Tube-only Experiment, a Plate Experiment, or an Experiment using imported files. This button is always visible and enabled except during acquisition.

**New Experiment from Template:** Opens the **New Experiment from Template** dialog (“New experiment from template dialog” on page 613), which enables you to create a new Experiment from an existing template with predefined settings. This button is always visible and enabled except during acquisition.

**Overlay Builder:** Opens the **Overlay builder** dialog (“Overlay Builder” on page 260), which enables you to select the samples and plots for inclusion in an Overlay plot. This button is only visible when the active experiment contains one or more samples and the experiment level workspace contains at least one plot. This button is not available during acquisition and it is disabled when a Compensation control is active.

### Templates group

**Templates** group enables you to save an active Experiment as a template.


**Save as Template:** Opens the **Save as Template** dialog (“Save As Template dialog” on page 747), which enables you to save the current Experiment as a template. This button is disabled on the Main Menu, Performance Test, and Instrument Configuration screens.

**Manage Templates:** Opens the **Manage Templates** dialog (“Manage templates dialog” on page 749), which enables you to edit, import, export, and delete templates. This button is disabled when acquisition is in progress.




## General group


**General** group enables you to customize the page layout and save an active Plate or Tube Experiment Workspace as a PDF.


 **Page Setup:** Opens the **Page Setup** dialog (“Page Setup dialog” on page 728), which enables you to customize the page layout (page size and orientation). This button is only available when Workspace, Reports, Results, Heat Map, or Overlay tab is active.

## Clipboard group

The **Clipboard** group enables you to cut, copy, and paste selected objects and text in the work area, and to Undo and Redo previous actions.

 **Paste:** Pastes the clipboard item into the current Workspace, Overlay gallery, Run Protocol, or Report. This option is only available when an item has previously been copied or cut into clipboard memory.

 **Cut:** Removes the selected item from the Workspace, Overlay gallery, or Report and retains it in memory until another item is cut or copied into memory. This option is only available when a Workspace, Overlay gallery, or Report item is selected.

 **Copy:** Retains the selected item in memory for pasting until another item is cut or copied in to memory. Workspace objects (including gates), Run Protocols, Overlays, Gallery plots, Report objects and Results can be copied into memory for pasting.

- Selecting **Copy** does not remove the item from the Workspace.
- When only a single statistics field, text field, or image is selected on the Workspace, the object is copied as a picture (enhanced metafile) to the Windows™ clipboard and as an HTML formatted object.

The object (stats field, text field, or image) can be pasted to an external application as a picture (enhanced metafile) or as HTML (CF\_HTML).


- When multiple workspace objects are selected, the selected objects are copied to the clipboard as an HTML formatted object and can be pasted to an external application as an HTML formatted object.


When objects are pasted as HTML formatted objects, the individual elements are pasted to the external application in the order they are selected.

- If plots are in the selection, the plots are copied as a picture (enhanced metafile) to the Windows™ clipboard and can be pasted as a picture to an external application.
- When Workspace objects are copied to the clipboard as HTML formatted objects, temporary files are created in:

C:\Users\\AppData\Local\Temp\AttuneNxTClipboard

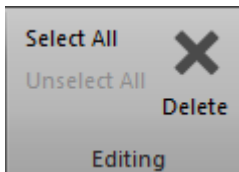
When the application closes, these temporary files are deleted.

 **Undo:** Negates the last undoable action in the order performed until the undo stack is empty. The **Undo** button is enabled only if there is an undoable action available. External functions including **Save**, **Import**, **Export**, or **Print** commands cannot be undone.

 **Redo:** Reverses the **Undo** action in the order performed until all available tasks have been carried out. This option is only available when an **Undo** action has been successfully performed.

## Editing group

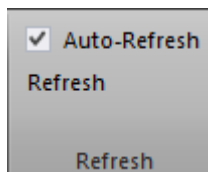
The **Editing** group contains shortcuts for selecting, unselecting, and deleting any of the objects in the work area.



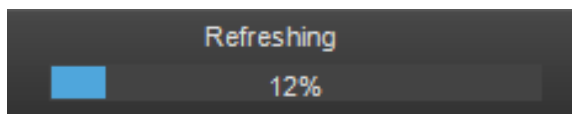
- **Select All:** Selects and highlights all objects on the current Workspace, Results, or Heat Map. This button is available if a Workspace, Results, or Heat Map is open and one or more items are present.
- **Unselect All:** Deselects all previously selected objects on the current Workspace. This button is available if a Workspace is open and one or more items are currently selected.
- **Delete:** Deletes the selected item. This button is available when a Workspace object or a Heat Map location is selected.

## Refresh group

The **Refresh** group contains options that enable the **Auto-Refresh** function to be turned off for inactive Workspaces, Heat Maps, and Results, therefore requiring them to be manually refreshed.



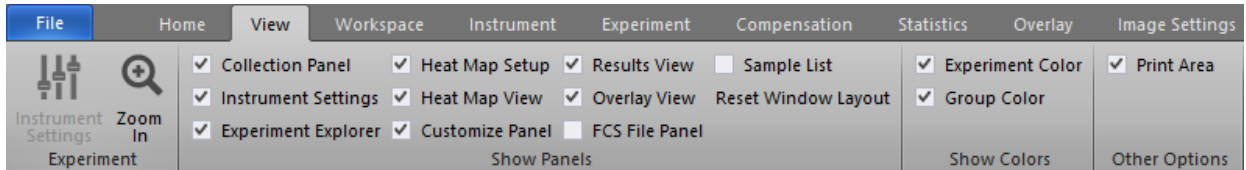
- **Auto-Refresh** checkbox: When **Auto-Refresh** is selected, all statistics in the Results tab are updated automatically in the background as the Workspace is modified. The Heat Map for all wells and tubes are also updated automatically.  
Deselecting **Auto-Refresh** turns off the automatic updates. However, printing and exporting statistics force an automatic refresh of all out-of-date items.  
The **Auto-Refresh** checkbox is enabled and checked by default.
- **Refresh:** Forces a refresh of all out-of-date items on the Results, Heat Map items, and Overlays. The **Refresh** button is always enabled.  
You can also access the **Refresh** function using the keyboard shortcut **F5**.



## View tab

**View** tab provides options to adjust the instrument settings and to control the overall look of the software. It is organized into four functional groups:

- **Experiment** (“Experiment group” on page 77)
- **Show Panels** (“Show Panels group” on page 78)
- **Show Colors** (“Show Colors” on page 79)
- **Other Options** (“Other Options group” on page 80)



## Experiment group

The **Experiment** group enables you to view and adjust Instrument Settings for an Experiment and to zoom in and out.

**Zoom In:** Displays all plots present in the **filmstrip area** of the **Experiment Workspace** tab (“Previews” on page 132) and enables you to zoom in (maximize) each plot. This is a toggle button (see “**Zoom Out**”) and is always visible; it is enabled when the **Experiment Workspace** is active and contains plots.

The plots are displayed in the **filmstrip area** in the order that they were created. When you click **Zoom In**, the plots are maximized according to these rules:

- If no plots are selected, the first plot is maximized.
- If one plot is selected, the selected plot is maximized.
- If more than one plot is selected, the first plot created is maximized.

**Zoom Out:** Closes the **Zoom In** function and restores the normal **Experiment Workspace**. This is a toggle button (see “**Zoom In**”) and is always visible; it is enabled when the **Experiment Workspace** is in the zoomed-in state.

---

**Note:** Alternatively, click the **Close** icon (X) to restore the normal **Experiment Workspace**.

---

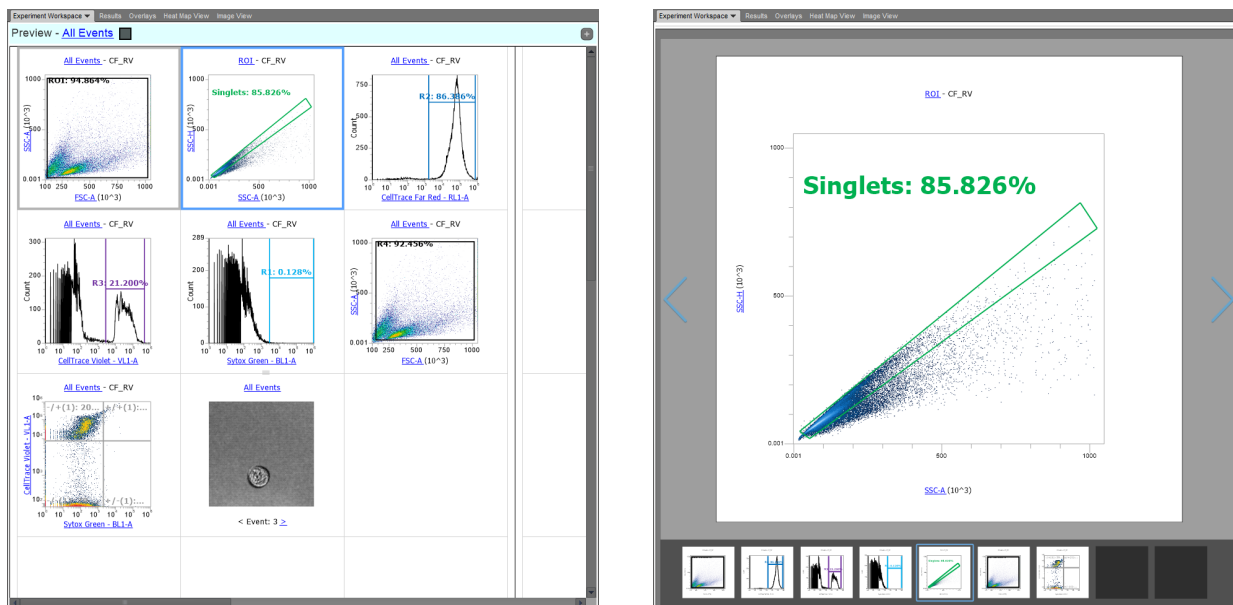
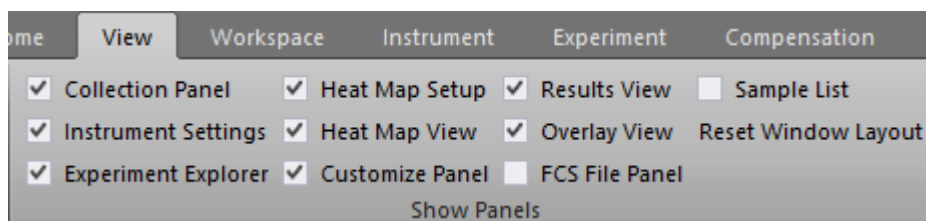


Figure 7 Experiment Workspace in the normal state (i.e., zoomed out-state) (left) and when Zoom In button is clicked (zoomed in-state) (right).

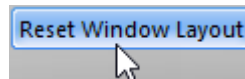
## Show Panels group

The **Show Panels** group enables you to open or close the panels and views in the **Main Application Workspace** (“Main application workspace” on page 56).



- **Collection Panel** (Chapter 12, “Collection panel”)
- **Instrument Settings** (Chapter 13, “Instrument settings panel”)
- **Experiment Explorer** (Chapter 11, “Experiment Explorer”)
- **Heat Map Setup** (Chapter 14, “Heat map setup panel”)
- **Heat Map View** (Chapter 6, “Heat Map View”)
- **Customize Panel** (Chapter 15, “Customize panel”)
- **Results View** (Chapter 8, “Results view”)
- **Overlay View** (Chapter 9, “Overlays”)
- **FCS File Panel** (Chapter 17, “FCS information panel”)
- **Sample List** (Chapter 7, “Sample List view”)
- By default, all options in the **Show Panels** group are always visible and enabled (except **FCS File Panel** and **Sample List**). For the default positions and display status of the panels and views, see “Docking behavior” under “Main Application Workspace” on “Docking locations” on page 58.

- To close a panel or view, deselect its checkbox.
- To open a panel or view, select its checkbox. Selecting a checkbox restores the panel or view to the state it was in when it was last closed.
- If a panel or view is hidden (“Auto Hide” on page 60), but not closed, its checkbox remains selected.
- By default, the **FCS File Panel** and **Sample List** are deselected, and all other panels and views are selected.
- Any changes you make for the **Show Panels** options persist for the next time you sign in.
- Click **Reset Windows™ Layout** to restore all panels and views to their default states.



---

**Note:** When hidden, the panel or view name appears as a tab on the side of the **Main Application Workspace** where the panel was docked.

---

## Show Colors

The **Show Colors** group contains the controls to show or hide the **Experiment** and **Group** colors in the **Heat Map View**.

- **Experiment Color:** When selected, shows the **Experiment color** for the **Samples** in the **Heat Map View**.  
When deselected, hides the **Experiment color** for the **Samples** in the **Heat Map View**.
- **Group Color:** When selected, shows the **Group color** for the **Samples** in the **Heat Map View**.  
When deselected, hides the **Group color** for the **Samples** in the **Heat Map View**.

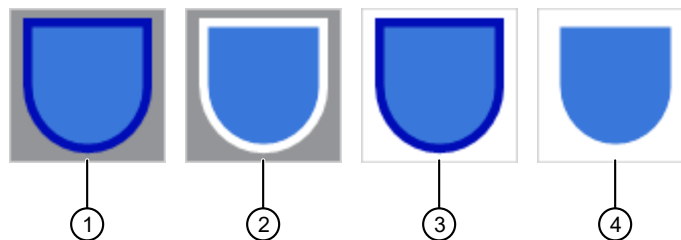


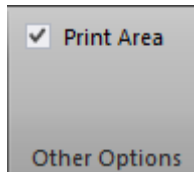
Figure 8 Effect of Show Colors controls on Samples in Heat Map View.

- ① **Experiment Color** and **Group Color** are both selected.
- ② **Experiment Color** is selected, **Group Color** is deselected.
- ③ **Experiment Color** is deselected, **Group Color** is selected.
- ④ **Experiment Color** and **Group Color** are both deselected.

## Other Options group

The **Other Options** group contains the display option for the **Print Area**.

The **Print Area** option is always visible and enabled.



### Print area

Enables you to show the Workspace view (Chapter 5, “Workspace view”) as it will be printed on a page.

- Select the **Print Area** option to show the page boundaries as gray lines drawn under an object.
- Deselect the **Print Area** option to restore the normal Workspace view.

## Workspace tab

**Workspace** tab provides tools for adding objects to a **Workspace** and aligning objects on the **Workspace**.

When using the Attune™ CytPix™ instruments or when the experiment supports imaging, the **Workspace** tab also contains the **Cell Image Analysis** tools.

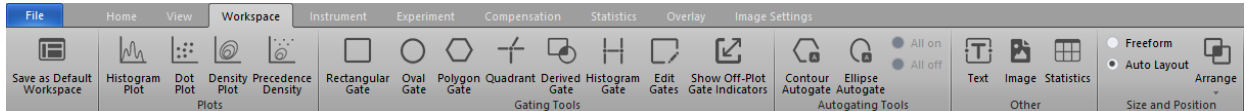


Figure 9 Workspace tab (Attune™ NxT Flow Cytometer)

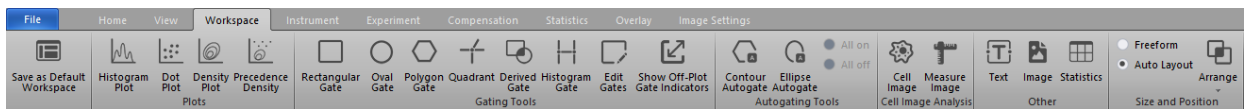


Figure 10 Workspace tab with Cell Image Analysis tools (Attune™ CytPix™ Flow Cytometer)

The **Workspace** tab is organized into these functional groups:

- **Save as Default Workspace** (“Save as Default Workspace” on page 81)
- **Plots** (“Plots group” on page 82)
- **Gating Tools** (“Gating Tools group” on page 83)
- **Autogating Tools** (“Autogating Tools group” on page 84)
- **Cell Image Analysis** (“Cell Image Analysis group” on page 85)
- **Other** (“Other group” on page 85)
- **Size and Position** (“Size and Position group” on page 86)

---


**Note:** **Cell Image Analysis** group is available only when an Attune™ CytPix™ Flow Cytometer is in use or when the Experiment supports imaging. This group is not visible for experiments created with the Attune™ NxT Flow Cytometer, which does not have imaging function.

---

**Note:** All options on the **Workspace** tab are inactivated during analysis of a **Compensation Control Sample**.

---

## Save as Default Workspace

 **Save as Default Workspace:** Saves the current Workspace as the default. Any new Experiments and Plates use this Workspace as the default Experiment Workspace.


The **Save as Default Workspace** option is inactivate when the Experiment Workspace (Chapter 5, “Workspace view”) is not in focus.


## Plots group


**Plots** group is used for creating plots of various types. **Plots** group controls are enabled when the Workspace view (Chapter 5, “Workspace view”) is in focus and at least two parameters are active, except for the Histogram plot control, which requires only one active parameter.




**Plots** group contains the following controls. For more information about plot types, see “Plots” under “Workspace view” (“Plots” on page 138).

 **Histogram Plot:** Adds a new single-parameter **Histogram plot** to the current Workspace. The Histogram plot displays the default parameter plotted against count (“Histogram plot” on page 139).

 **Dot Plot:** Adds a new **Dot plot** to the current Workspace. The Dot plot displays two-parameter data where each axis represents the signal intensity of one parameter. Different colors are used to represent the parent gate of events that fall within bins on the plot (“Dot plot” on page 140).

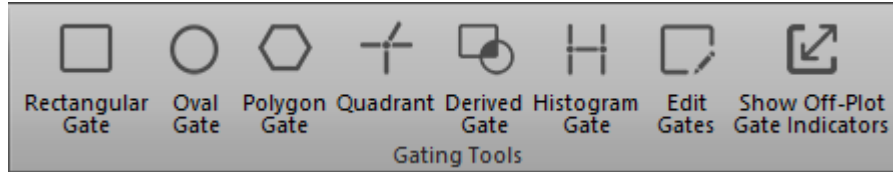
 **Density Plot:** Adds a new **Density plot** to the current Workspace. The Density plot displays two-parameter data where the colors represent the collection of events with the same intensity and each axis represents the signal intensity of one parameter (“Density plot” on page 141).

 **Precedence Density Plot:** Adds a new **Precedence Density plot** to the current Workspace. The Precedence Density plot is a combination of Dot and Density plots, where a gradient is used to indicate the number of events within each plot bin, and color is used to indicate the parent gate of events present (“Precedence density plot” on page 142).



## Gating Tools group


**Gating Tools** enable you to isolate a region in a selected plot for analysis. **Gating Tools** are available when the Workspace view (Chapter 5, “Workspace view”) is displayed, unless otherwise specified.



- Dual-parameter gate types are active when the current Workspace contains at least one dual-parameter plot (i.e., Dot plot, Density plot, Precedence Density plot).
- Single-parameter gate types are active when the Workspace contains at least one Histogram plot.
- A maximum of 128 gates can be created on plots on a Workspace, except Quadrant gates, of which at least 1,000 can be created. The Gating Tools controls are inactive when the Workspace reaches the maximum number of gates allowed.

The **Gating Tools** group contains the controls listed below. For more information about gates, including instructions for creating and deleting gates, see “Gates” under “Workspace view” (“Adjust plots” on page 143).

 **Rectangular Gate:** Enables you to draw a **Rectangular gate** on a dual-parameter plot as described in “Rectangular Gate” on page 149.


 **Oval Gate:** Enables you to draw an **Oval gate** on a dual-parameter plot as described in “Oval Gate” on page 150.

 **Polygon Gate:** Enables you to draw a **Polygon gate** on a dual-parameter plot as described in “Polygon Gate” on page 151.

 **Quadrant Gate:** Enables you to draw a **Quadrant gate** on a dual-parameter plot as described in “Quadrant Gate” on page 152.

 **Derived Gate:** Displays the **Derived Gate dialog**, which enables you to create gates based on Boolean operators (i.e., **Derived gates**) as described in “Derived gate” on page 163. **Derived Gate** control is active when the Workspace view is displayed and the current Workspace has at least one gate present.

 **Histogram Gate:** Enables you to draw a **Histogram gate** on a Histogram plot as described in “Histogram Gate” on page 155.

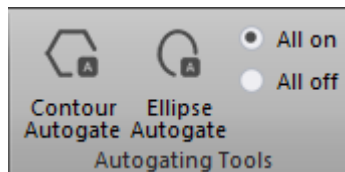
 **Edit Gates:** Displays the **Edit Gates dialog**, which enables the editing of gate color, parent gate, gate math expression, the z-order in which gates are painted as described in “Edit gates dialog” on page 742. **Edit Gates** control is available when the Workspace view is displayed and the current Workspace has at least one gate present.

 **Show Off-Plot Gate Indicators:** Displays **Off-Plot Gate Indicators** in plots that include gates that exist outside of the plot limits as described in “Off-plot gates” on page 164.

## Autogating Tools group


**Autogating Tools** enable you to create auto region bounding rectangles in a selected plot, which determine the autogate target area when identifying populations based on autogate settings. If the criteria defined in autogate settings are met, the target area becomes autogated (see “Autogate gating rules” on page 162). Autogates function in the same way as regular gates to isolate a region for analysis.

**Autogating Tools** are available when the Workspace view (Chapter 5, “Workspace view”) is displayed, unless otherwise specified.

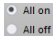


- **Autogating Tools** are enabled when the current Workspace contains at least one dual-parameter plot (i.e., Dot plot, Density plot, Precedence Density plot). If the Workspace does not contain any plots, then all buttons in the **Autogating Tools** group are disabled.
- A maximum of 128 gates, including autogates, can be created on plots on a Workspace, except Quadrant gates, of which at least 1,000 can be created. The **Autogating Tools** controls are disabled when the Workspace reaches the maximum number of gates allowed.

The Autogating tools group contains these controls:

 **Contour Autogate:** Enables you to create a **Contour auto region** on a dual-parameter plot as described on “Contour Autogate” on page 156.

 **Ellipse Autogate:** Enables you to create an **Ellipse auto region** on a dual-parameter plot as described on “Ellipse autogate” on page 159.

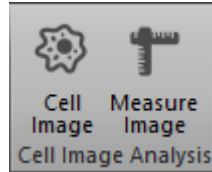
 **All on and All off:** The **All on** button turns all autogated and unautogated regions in the currently active Workspace ON. The default state for all autogates is to be in an ON state with the **All on** button selected.

The **All off** button turns all autogated and unautogated regions in the currently active Workspace OFF.

The **All on** and **All off** button states persist with the Workspace.


## Cell Image Analysis group


**Cell Image Analysis** group enables you to add a new **Cell Image Container** to the Workspace and to perform size measurements on the image displayed in the image container.



The **Cell Image Analysis** tools are available only when using an Attune™ CytPix™ Flow Cytometer or when the Experiment supports imaging. This group is not visible for experiments that were created using an Attune™ NxT Flow Cytometer, which does not have imaging capability.

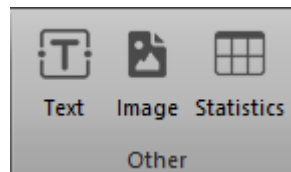
The **Cell Image Analysis** group contains these controls:

 **Cell Image:** Adds a new **Cell Image Container** to the current Workspace, which enables you to add a cell image specific to the Sample (“Cell image container” on page 181). Cell image containers act like plots and can be batch printed with the Workspace (cell images cannot be printed from the Image view Image Gallery).

 **Measure Image:** Lets you to draw a measurement tool on a cell image container to measure the area of the cell in the image (“Measure Image tool” on page 184). This control is only enabled if the Workspace contains a cell image container.

## Other group


**Other** group is used to insert text and images into the Workspace, and to show the statistics that correspond to a selected plot in the Workspace.



The **Other** group contains the following controls, which are available when the Workspace view is in focus.

 **Text:** Adds a new **Text box** to the current Workspace.

- The **Text box** displays the default text “New Text Box”, which can be edited in place without opening a separate dialog box.
- The style and border of the **Text box** can be customized as described in “Customize Text Box Options” on “Customize text box options” on page 444.

 **Image:** Opens a standard Windows™ file selection dialog for selecting an image file to add to the Workspace.

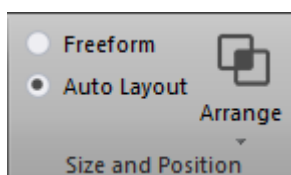
- The image file cannot be larger than one page and is scaled to fit on the page preserving the aspect ratio, if needed.
- The image types available for insertion include JPG, GIF, BMP, PNG, TIF, and Windows™ enhanced metafile.

 **Statistics:** Adds a new **Statistics box** to the current Workspace.

- If no plot is selected, clicking on the **Statistics** button inserts a **Workspace statistics table** in the default Workspace location.
- If one or more plots are selected, then one **Plot Statistics box** is added to each selected plot immediately below the relevant plots as described in “Workspace statistics” on page 175.

## Size and Position group

The **Size and Position** group provides tools that assist in creation and customization of Workspaces, and contains the buttons listed below.



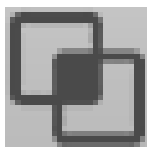
### Layout

**Layout** controls consist of two radio buttons, which control the **View mode** of the Workspace layout as described in “Workspace view modes” on “Workspace view modes” on page 126.

- **Auto Layout:** In this mode, objects are placed into grid slots. The number of available slots on the Workspace is defined by the **Grid Size** options (“Grid options” on page 87). **Auto Layout** is the default mode.
- **Freeform:** In this mode, objects can be freely resized and moved anywhere on the Workspace. There is no automatic arrangement or alignment of objects on the Workspace.

### Arrange

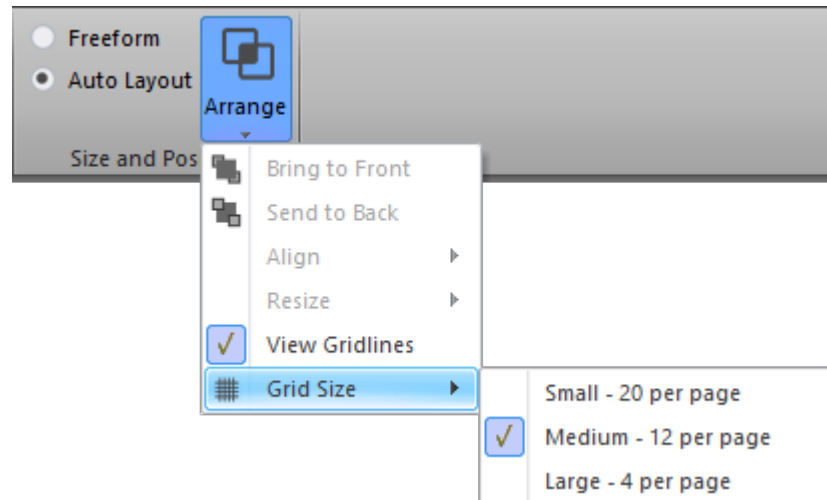
**Arrange** opens the **Arrange Tools** dropdown menu, which contains the tools to arrange, align, and size multiple Workspace objects.



- The **Arrange Tools** commands are available only when multiple items are selected on the Workspace, unless otherwise stated.
- The **View Gridlines** and **Grid Size** options are only enabled in the **Auto Layout** mode; all other options are enabled only in the **Freeform** mode.

## Grid options

**Grid options** are enabled in the **Arrange Tools** dropdown menu only when the **Auto Layout** mode is selected using the **Layout** radio buttons (“Grid options” on page 87).



**View Gridlines:** Selecting this option displays the lines that form the grid onto which the Workspace objects are snapped. By default, the gridlines are displayed.

**Grid Size:** Opens the **Grid Size** dropdown menu, which contains the options to define the number of grid slots available per Workspace page. A check mark appears next to the currently selected size.

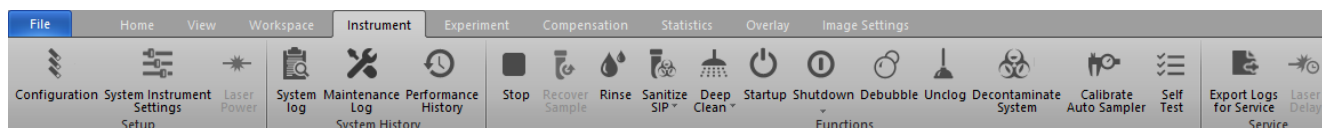
The default selection is 12 grid slots per page.

You can select a larger grid size allowing only 4 items per page, or a smaller grid size allowing up to 20 items per page at the default sizes.

## Instrument tab

**Instrument** tab enables you to view and export instrument performance and maintenance data, to update instrument configuration, to view and change system instrument settings, and to perform instrument maintenance tasks. It is organized into five functional groups:

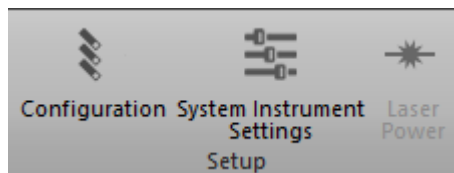
- **Setup** (“Setup group” on page 88)
- **System History** (“System History group” on page 89)
- **Functions** (“Functions group” on page 90)
- **Service** (“Service group” on page 93)
- **Automation** (“Automation group” on page 94)



**Note:** Most buttons in the **Instrument** tab are inactive in **Automation** mode (“Automation group” on page 94), with a few becoming active when Plate Acquisition is paused in Automation mode.

### Setup group

The **Setup** group enables you to view instrument data and manage configuration files.



**Configuration:** Opens the **Configuration module** (“Filter configuration (FC) module” on page 512), which enables you to manage instrument configuration files and filters. This button is always visible and enabled except during acquisition.

**System Instrument Settings:** Opens the **System Instrument Settings** dialog (“System instrument settings” on page 401), which enables you to modify **Area scaling factor (ASF)** and **Window extension** settings.

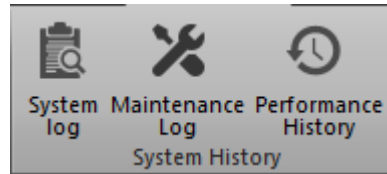
This button is visible only to users that have privileges to modify **System Instrument Settings**.

Access to **System Instrument Settings** is disabled when the instrument is acquiring a sample or a plate, running Baseline/PT, while paused, and when set to automation mode.

**Laser Power:** Enables users with Service accounts to turn the lasers on and off. This button is available only for Service accounts (i.e., Thermo Fisher Scientific service team); it is invisible to all other account types.


## System History group

The **System History** group enables you to view the **System Log**, **Maintenance Log**, and **Performance History** reports.



 **System Log:** Opens the **General** tab of the **System Log dialog** (“System Log dialog” on page 793), which enables you to view system transactions.

- The information displayed in the **System Log** depends on the type of account that you have: service, administrator, system administrator, or general user account.
- The **System Log** button is always visible and active.

 **Maintenance Log:** Opens the **Maintenance** tab of the **System Log dialog** (“Maintenance tab” on page 800), which lists all the maintenance functions that were performed on the instrument.


- The **Maintenance** tab of the **System Log dialog** contains a system generated record of all the instrument maintenance tasks performed by all users using the **Functions** group of the **Instrument** ribbon tab. The record also includes the start and stop dates and times of these tasks.
- The **Maintenance Log** also enables you to log other maintenance actions that are not listed on the **Instrument** ribbon tab, such as filter replacements.
- The **Maintenance Log** can be sorted by date, type, state, and user, and can be exported locally or printed to keep record of the instrument history.
- The **Maintenance Log** button is always visible and active.

---

**Note:** In the SAE mode, you need to have **Attune Software > Other > View system logs** permission selected to view the **Maintenance** tab and to add custom maintenance tasks (see “SAE account permissions” on page 887).

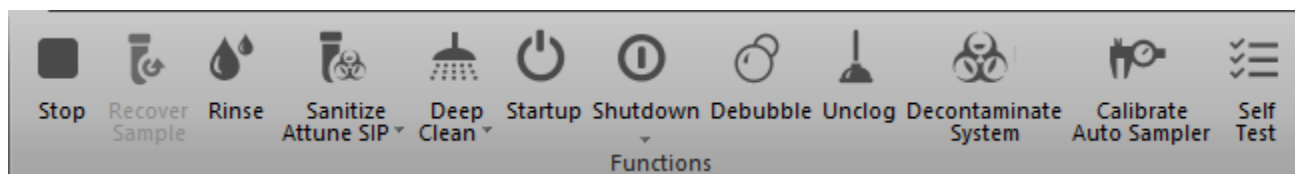
Both the **Security Configuration** and the **Audit History** permissions must be selected in the **SAE Administrator Console** (see “SAE account permissions” on page 887) to see the **User** column in the **Maintenance Viewer**, and to view which user has performed which instrument function.

---

 **Performance History:** Opens the **Performance History Report** (“Performance History Report” on page 557), which enables you to view the pass/fail status of all Performance Tests. This button is always visible and active except during acquisition.

## Functions group

The **Functions** group enables you to start the instrument functions.



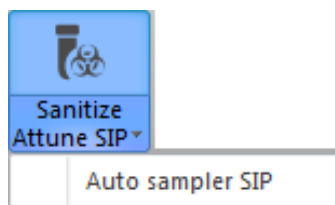
**Stop**: Stops any acquisition or instrument function, except the **Calibrate**, **Self Test**, and **Rinse** instrument functions. This button is always visible and active.

- If you click **Stop** during the following instrument functions, you need to run **Startup** before running the instrument again:
  - **Startup** (Startup on page 91)
  - **Unclog** (Unclog on page 91)
  - **Shutdown** (Shutdown on page 91)
  - **Deep Clean** (Deep Clean on page 91)
  - **Decontaminate System** (Decontaminate System on page 91)
- If you click **Stop** during the **Rinse** function, the instrument ignores the **Stop** command.
- If you click **Stop** when running a Plate Sample, the instrument stops the acquisition and aborts the Plate.
- If you click **Stop** during all other instrument functions, the instrument forces a **Rinse** before performing any other operation.
- Clicking **Stop** when the instrument is in the Automation mode toggles the Automation mode to 'OFF'.

**Recover Sample**: Recovers the unused sample from the sample loop. This button is active when the instrument status is idle (that is, it is not currently acquiring) and there is sufficient sample in the sample loop or sufficient preloaded sample to recover. When the button is pressed, the remaining sample is returned into a tube (from the sample loop) or back into the sample well (preloaded sample in a plate).


**Rinse**: Opens the **Rinse** dialog, which enables you to rinse the Sample lines. This button is always visible; it is active when the instrument status is idle and not currently acquiring a plate.

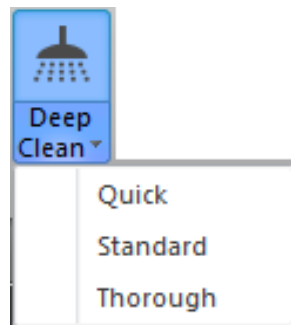
**Sanitize Attune™ SIP** : Opens the **Sanitize** dialog, which enables you to sanitize the instrument SIP (sample injection port) and sample lines, or the Auto Sampler SIP and Sample lines. This button is always visible; it is enabled when the instrument status is idle and not actively acquiring.




This is a split button: Click **Sanitize Attune™ SIP** or select **Auto Sampler SIP** from the dropdown list.




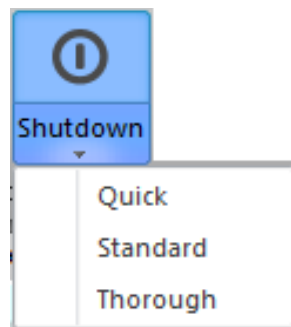
 **Deep Clean:** Opens the **Deep Clean** dialog, which enables you to thoroughly wash the system sample lines and flow cell. This button is always visible; it is active when the instrument status is idle and not currently acquiring. It is inactive when a Plate run is paused.




This is a split button: Click **Deep Clean** to perform a standard clean cycle, or select **Quick**, **Standard**, or **Thorough** from the dropdown list.


 **Startup:** Opens the **Startup** dialog, which enables you to prime the system fluidics. This button is always visible; it is active when the instrument status is idle and not currently acquiring, or when the instrument is in the “sleep” state. It is inactive when a Plate run is paused.


 **Shutdown:** Opens the **Shutdown** dialog, which enables you to sanitize and shut down the instrument. This button is always visible; it is active when the instrument status is idle and not currently acquiring. It is inactive when a Plate run is paused.



This is a split button: Click **Shutdown** to perform a standard shutdown cycle, or select **Quick**, **Standard**, or **Thorough** from the dropdown list.


 **De-bubble:** Opens the **De-bubble** dialog, which enables you to clear the sample lines and flow cell of bubbles. This button is always visible; it is active when the instrument status is idle.

 **Unclog:** Opens the **Unclog** dialog, which enables you to unclog the Sample lines. This button is always visible; it is active when the instrument status is idle and not currently acquiring.

 **Decontaminate System:** Opens the **Decontaminate** dialog, which enables you to decontaminate all system fluid lines. This button is always visible; it is active when the instrument status is idle and not currently acquiring. It is inactive when a Plate run is paused.

If an Attune™ Auto Sampler is connected to the system and powered on, **Decontaminate System** initiates a decontamination cycle for the cytometer and the Auto Sampler.


For detailed instructions about instrument decontamination, see the *Attune™ Flow Cytometry Maintenance and Troubleshooting Guide* (Pub. No. 100024234), available for download at [thermofisher.com](http://thermofisher.com).

 **Calibrate Auto Sampler:** Opens the **Auto Sampler Calibration** dialog, which enables you to calibrate the Attune™ Auto Sampler. **Calibrate Auto Sampler** button is visible only when there is an auto sampler connected to the instrument. The button is active when the instrument status is idle and not currently acquiring; it is inactive when a Plate run is paused.

---

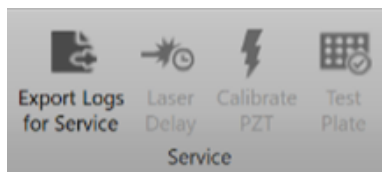
**Note:** Auto Sampler calibration automatically occurs every 30 days in Attune™ Cytometric Software versions 2.5 or greater.

---

 **Self Test (System Test):** Opens the **System Test** dialog, which lets you run system diagnostics. This button is visible only to Service accounts; it is active when the instrument status is idle and not currently acquiring. It is inactive when a Plate run is paused.


## Service group


The **Service** group contains functions for servicing the instrument.



**Laser Delay** and **Calibrate PZT** functions are available only to users with Service accounts (i.e., Thermo Fisher Scientific service team).

**Export Logs for Service** function is available to all users.

 **Export Logs for Service:** Opens the **Export Logs for Service** dialog (“Export Logs for Service dialog” on page 772), which enables you to export selected logs to a single zip file based on the specified date range. This button is enabled, if the instrument status is idle or when a Plate run is paused. It is disabled, when moving between wells.

 **Test Plate:** Opens the **Test Plate** dialog (“Test Plate dialog” on page 778), which enables you to validate a plate to ensure that the autosampler probe position is in the correct location in all four corners of the plate.


The **Test Plate** button is only visible if the system is connected to or set to a CytKick™ Autosampler and the autosampler is powered on.

The **Test Plate** button is enabled if the instrument status is idle and an autosampler is connected and powered on. This function is disabled if the instrument is paused or set to automation mode.


## Automation group

The **Automation** group lets you to put the application under the control of external software to enable the automated loading and running of plates. This group is only visible if a TCP/IP Communication port has been specified in the command line (i.e., `attune.exe /p:8225`) or in the **Configuration Options** dialog (see “Automation communication settings” on page 697).



 **Connect Robot:** Sets the software in **Automation mode**, which puts the application under the control of external software. This button is only enabled if there is an Auto Sampler present and the instrument status is idle.

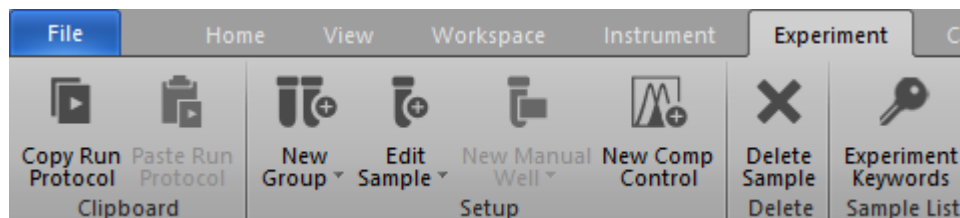
- When the **Connect Robot** button is toggled ON, the software is set to **Automation mode** and the button is highlighted blue.
- When in **Automation mode**:
  - All fluid functions are disabled, the Experiment cannot be changed, and navigating away from the Workspace view (i.e., Attune™ desktop) is disabled:
  - All **File Menu** commands except for **Exit** are disabled.
  - The **Instrument** tab **Setup** group buttons are all disabled unless a plate run is paused.
- If the instrument is disconnected from the external device, the toggle state is automatically turned OFF and the **Connect Robot** button is disabled.
- When the **Connect Robot** button is toggled OFF, the software accepts normal user inputs.

 **Map Plates:** Opens the **Plate Mapping** dialog (“Plate Mapping dialog” on page 756), which enables you to map defined templates to barcodes to be used when in **Automation mode**.

## Experiment tab

**Experiment** tab provides tools to define the Experimental layout for tubes and plates. The **Experiment** tab is visible only when the **Heat Map view** (Chapter 6, “Heat Map View”) is active.

Any changes that you make to the Experiment layout in the **Experiment** tab are reflected in real-time in the **Experiment Explorer** (Chapter 11, “Experiment Explorer”).



## Experiment limits


Plate-based Experiments can contain Samples in Wells and/or Tubes. Unless otherwise specified, in this section Sample is used to mean Samples in Tubes and in Qells.

All Samples on a Plate Experiment must total  $\leq 400$  Tells and tubes. This limits the number of Tube Samples per Plate to:


- $\leq 304$  Tube Samples for 96-well plates
- $\leq 16$  Tube Samples for 384-well plates

Changes made to the Experiment in the **Heat Map view** are automatically reflected in the **Experiment Explorer**.

## Experiment tab controls

 **Copy Run Protocol:** Copies the **Run Protocol** for the selected samples to the clipboard. This button is always visible, but it is enabled only when samples are selected.

- If you select multiple samples that have different Run Protocols, the software copies all protocols and maintains the relative positions of the selected samples.

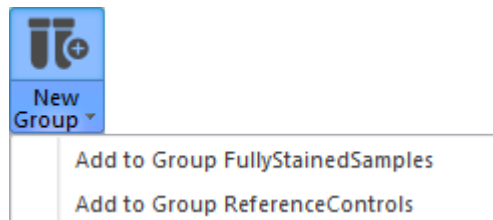
 **Paste Run Protocol:** Pastes **Run Protocol** from the clipboard into the selected samples. This button is always visible, but it is enabled only when one or more Run Protocols are copied to the clipboard and samples are selected.

- If you copy only one Run Protocol to the clipboard, the software pastes that Run Protocol into all selected samples.
- If you copy multiple Run Protocols to the clipboard, this button is active only when the same array of samples is selected. The software pastes each unique Run Protocol into the selected sample of the same position in the array.
- If the Run Protocol is copied from and pasted into plates of different types, the software checks all entries for validity following the rules set in “Run Protocol Options” (“Record dialog for manual well” on page 367).

For example, if the pasted well volume is greater than allowed in the selected well, the software defaults to the maximum volume allowed.

**New Group:** Adds the selected Samples to a new or existing Group. This button is always visible, but it is enabled only when samples are selected. The selected samples can be unprogrammed or marked as part of the Experiment.

The **New Group** button is a split button that includes a dropdown menu option.



- To create a new Group and add the selected Samples to it, click the **main portion** of the **New Group** button.
- To add the selected Samples to an existing Group, select the **Group** from the **New Group dropdown** menu.

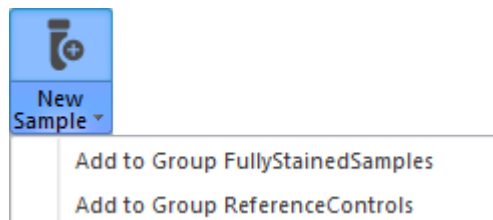
---

**Note:** Each **Experiment** that has assigned **Samples** consists of one or more **Groups**. Each **Group** can contain up to 400 **Samples**, and **Sample** can be derived from a **Well** in a **Plate** or from a **Tube**.

---

**New Sample:** Marks selected blank tubes or wells as Samples and adds these new Samples to a Group. This button is always visible, but it is enabled only when blank tubes and/or wells are selected.

The **New Sample** button is a split button that includes a dropdown menu option.



- To add the new Samples to the most recently created Group, click the **main portion** of the **New Sample** button. If no groups exist, the software creates a new Group.
- To add the new Samples to an existing Group, select the **Group** from the **New Sample dropdown** menu.

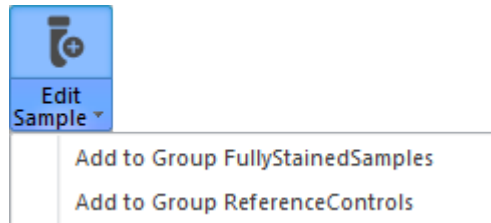
---

**Note:** The **New Sample** button is a toggle button, which changes to **Edit Sample** when one or more Wells and/or Tubes already marked as Samples are selected (see “Edit Sample” on page 96).


---

**Edit Sample:** Moves selected Samples to a different Group. This toggle button (see “**New Sample**” above) is always visible, but enabled only when Samples are selected.

The **Edit Sample** button is a split button that includes a dropdown menu option.

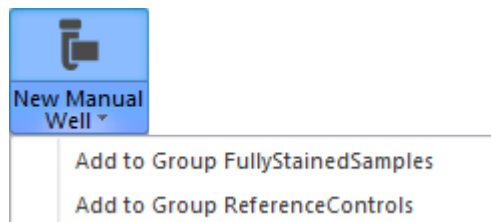


- To move the selected Samples to a different Group, select the **Group** from the **Edit Sample dropdown** menu.
- If the selection also includes new Samples, the button remains **Edit Sample**. New Samples are added to the same Group selected for the existing Samples.

 **New Manual Well**: Creates a new Sample in the selected wells marked as Manual Wells in the most recently created Group. If no Groups exist, a new Group is created and the Manual Well samples are added to this group.

The **New Manual Well** button is always visible, but it is enabled only when one or more wells are selected. The button is disabled for Tube-based Experiments or if Tube locations are selected on the Heat Map in a Plate-based Experiment.

The **New Manual Well** button is a split button that includes a dropdown menu option.



- The software marks Manual wells with an **MW** icon  in the top left corner of the well.



- The tooltip for the Manual well button reads “Defines a well that has both ‘Run’ and ‘Record’ modes.”

 **New Comp Control**: Opens the **Compensation Setup** dialog (“Compensation setup dialog” on page 581).

- **Compensation Setup** dialog lets you select the compensation source from wells or tubes. If blank wells are selected before clicking **New Comp Control**, the compensation dialog selects wells as the compensation source and the selected wells are prepopulated with the Compensation controls. By default, all possible Compensation controls are selected. Note that each parameter selected for compensation needs to have an associated well location before the dialog is closed.

- Compensation wells are indicated in the software by the **CS** icon in the top left corner of the well. The well is also labeled with the fluorescence channel or the unstained control that is mapped to the compensation well.



**✕ Delete Sample:** Deletes the selected Samples, which returns the Samples to the undefined state and removes all associated sample data (Workspaces, Run Protocols, and FCS files). This button is always visible, but it is enabled only when Samples marked as part of an Experiment are selected.

**🔑 Experiment Keywords:** Opens the **Experiment Keywords** dialog (“Experiment Keywords dialog” on page 770), which enables you to add, remove, or edit experiment keywords. This button is only enabled when an Experiment is active, the Workspace is in view, and the system is not acquiring, paused, or in the Automation mode.

## Hot keys

Hot keys can be used alone or in combination to copy and paste Run Protocols and delete selected Samples.

- **Ctrl+C** combination on the keyboard copies the selected Run Protocols to the clipboard as described in “Copy Run Protocol” (“Experiment tab controls” on page 95).
- **Ctrl+V** combination on the keyboard pastes the Run Protocol in the clipboard to the selected Sample as described in “Paste Run Protocol” (“Experiment tab controls” on page 95).
- **Delete** key on the keyboard deletes selected Samples as described in “Delete Sample” (“Experiment tab controls” on page 95).



## Compensation tab

**Compensation** tab contains controls for setting compensation, selecting the compensation source, and changing the view of Compensation samples.

The contents of the **Compensation** ribbon tab vary depending on whether a **Sample Workspace** or a **Compensation Workspace** is displayed.

Depending on the context, the **Compensation** tab contains the following functional groups:

- **Setup** (“Setup group” on page 99)
- **Apply** (“Apply group” on page 100)
- **Adjustment** (“Adjustment group” on page 101)

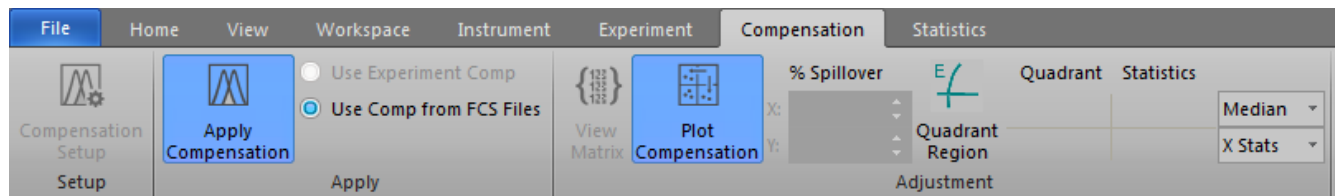


Figure 11 Compensation tab fully expanded on a Compensation Workspace.

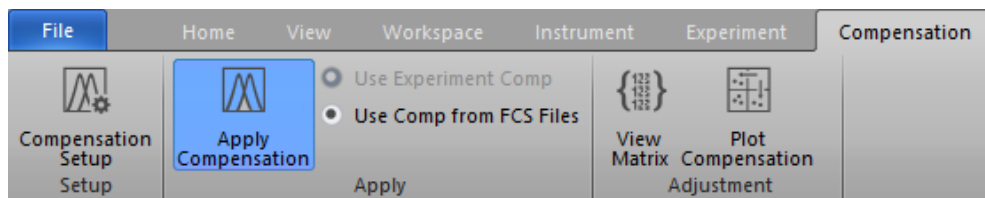
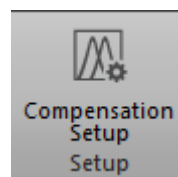


Figure 12 Compensation tab fully expanded on a Sample Workspace.

**Note:** For more information about the compensation function of the Attune™ Cytometric Software, see Chapter 21, “Compensation”.

## Setup group

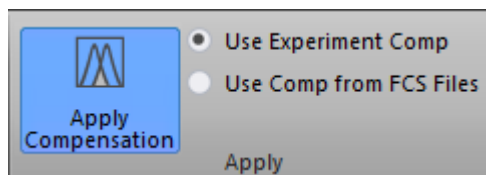
**Setup** group contains the **Compensation Setup** button. When the Compensation Workspace is active, more controls are visible that enable you to select **Histogram View**. The **Setup** group is always visible.




 **Compensation Setup:** Opens the **Compensation Setup dialog** as described on “Compensation setup dialog” on page 581. The **Compensation Setup** button is always visible, but it is enabled only when a Plate or Tube Experiment is active.

## Apply group

**Apply** group includes controls to turn compensation on and off for the selected Workspace and to select between Experiment Compensation or compensation loaded from the FCS headers. The **Apply** group is only visible on a Sample Workspace.



 **Apply Compensation:** When selected, all data shown on the Workspace and used for calculating statistics is compensated.

When deselected, all data shown on the Workspace and used for calculating statistics is the raw uncompensated data.

- The **Apply Compensation** button is **ON** by default, if compensation is available. The button state persists on a per Experiment basis.
- If Experiment-level compensation or FCS file compensation do not exist, **Apply Compensation** opens the **Compensation Setup dialog** (“Compensation setup dialog” on page 581).

---

**Note:** The **Apply Compensation** button is a state button; it remains **ON** until turned **OFF**. When active, the selected compensation option (**Use Experiment Comp** or **Use Comp from FCS Files**) is applied. If an alternative compensation option is selected, the settings are updated automatically.

---

**Use Experiment Comp** and **Use Comp from FCS Files** options are only visible and enabled when a Sample is active. These options are disabled on the Compensation Workspace and when compensation is not available from either the Experiment or the selected FCS file.

The button state persists on a per Experiment basis except for Samples where the selected compensation is invalid.

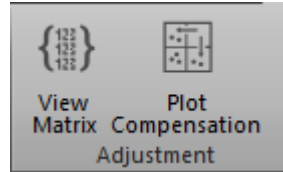
**Use Experiment Comp:** Uses the **Compensation Matrix** that is defined using Compensation controls or an imported compensation settings file. This is the default selection when Experiment-level compensation exists.

**Use Comp from FCS Files:** Uses the **Compensation Matrix** values stored in the FCS header. This is the default when Experiment compensation does not exist

- If an FCS file does not contain a **Compensation Matrix** and the **Use Comp from FCS Files** option is selected, then the **Use Experiment Comp** option is used for that Sample, if available.
- If an Experiment-level compensation is incompatible with an FCS file in the Experiment (due to importing an FCS file), then **Use Comp from FCS Files** option is used for that Sample, if available.
- In these cases, the invalid option is disabled. If the selected option is disparate to the Compensation used for a selected Sample (due to the Sample-level override), the buttons show the indeterminate state.

## Adjustment group

**Adjustment** group includes controls to view the **Compensation Matrix** and to enable the **Plot Compensation** function.



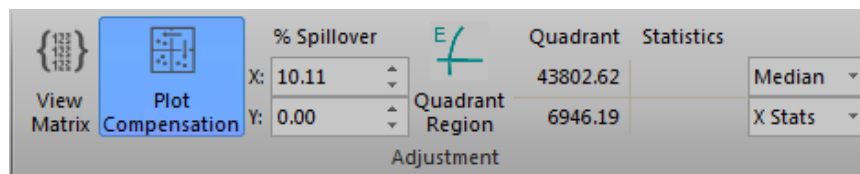
**View Matrix:** Opens the **Matrix dialog** (“Matrix dialog” on page 599), which shows the active compensation associated with the current Experiment or file. The **View Matrix** button is always visible, but it is only enabled when a Plate or Tube Experiment is active and either Experiment-level compensation settings exist or there is an active Sample containing an FCS file.

**Plot Compensation:** Enables **On-Plot Compensation Adjustment** (“On-Plot compensation adjustment” on page 602) and controls the visibility of the **Compensation Adjustment Tools** (below).

- The **Plot Compensation** button is deselected (**OFF**) by default.
- The button is enabled and visible only when **Apply Compensation** is selected. It is not visible in the Compensation Workspace, unless the **View Results** button is selected.
- Any changes made to compensation persist for both the Experiment-level compensation and the FCS compensation.
- The **Plot Compensation** button is a state button; it remains **ON** until it is turned **OFF** or if the selection of Workspace or the Sample changes.

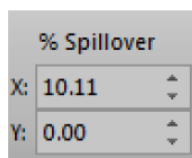
## Compensation Adjustment tools

- When **Plot Compensation** is turned **ON**, a collection of compensation adjustment tools are displayed on the **Compensation** tab.



- The **Compensation Adjustment tools** enable the manual adjustment of a selected plot, the ability to add a **Quadrant gate** to a plot, and the show **Quadrant statistics** as described below.
- When **Plot Compensation** is turned **OFF**, these controls are not visible.


**% Spillover Adjustment:** These tools consist of two controls that enable the adjustment of the compensation or the spillover matrix.



- These controls are only enabled when a dual-parameter plot with two fluorescence parameters is selected in a Sample Workspace or on the Compensation Workspace when **View Results** is turned **ON**.
- The top control corresponds to the X-axis and the bottom control corresponds to the Y-axis. The **up** and **down arrows** of the control or the keyboard move the spillover matrix value by increments of 1 with each click or continuously by maintaining the click.
- Pressing the **Shift** key when clicking on the **arrows** or when using the keyboard **up/down arrow** buttons moves the value by increments of 0.1.
- The text field accepts numbers from 0 to 100 and displays 2 decimal places.

 **Quadrant Gate:** Enables the insertion of a **Quadrant gate** on a dual-parameter plot as described on “Gating Tools group” on page 83.

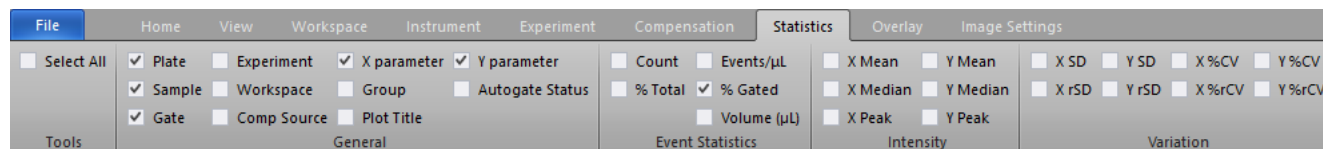
**Quadrant Statistics:** Enables the selection of the mean or the median values and the Y-axis or X-axis statistics.

	Quadrant	Statistics
	43802.62	Median ▾
Quadrant Region	6946.19	X Stats ▾

- By default, the **Median** and the **X Stats** (X-axis statistics) are selected.
- The statistics values are displayed within a representation of a quadrant gate. When selected, the X-axis statistics are shown in the lower left and lower right quadrants, and the Y-axis statistics are shown in the lower left and upper left quadrants.
- The values shown are based on the quadrant set on the currently selected plot.
- For the statistics to be displayed, a single dual-parameter plot containing only fluorescence parameters (not scatter) must be selected, and the plot must show a quadrant. In addition, an FCS file that can be compensated must be present. If these requirements are not met, the control does not show any statistics.
- If a statistics value cannot be calculated, the respective quadrant shows **N/A** (i.e., not applicable).

## Statistics tab

**Statistics** ribbon tab enables you to select which statistic is displayed on the Plots, Workspaces, and the Results view. The statistics are grouped in the ribbon according to their statistical type (**General**, **Event Statistics**, **Intensity**, and **Variation**).



The Statistics tab contains the following functional groups:

- **Results** (“Results group” on page 105)
- **Tools** (“Tools group” on page 106)
- **General** (“General group” on page 107)
- **Event Statistics** (“Event Statistics group” on page 108)
- **Intensity** (“Intensity group” on page 108)
- **Variation** (“Variation group” on page 109)

---

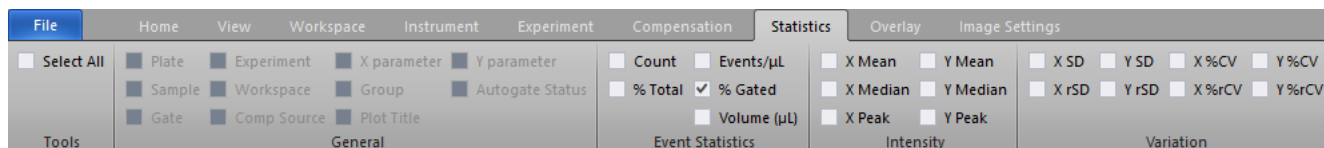
**Note:** The **Statistics** tab behaves differently (as explained below) depending on which Application tab is in focus or which object in a Workspace is selected. The entire **Statistics** tab is disabled when the **Heat Map View** is in focus (Chapter 6, “Heat Map View”).

---

## Statistics tab behavior

### Behavior in Workspace view

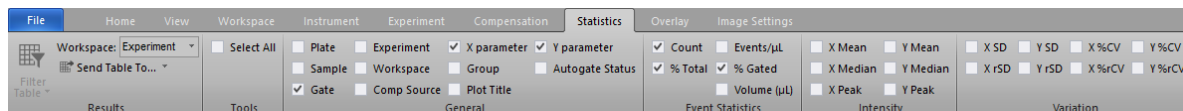
When the **Workspace** view (Chapter 5, “Workspace view”) is in focus and a statistics table is **not** selected, only a single statistic can be selected at a time. This selection applies to the statistic displayed on all plots. In the **Workspace** view, the **Results** group (“Results group” on page 105) is not visible and the **General** group (“General group” on page 107) is not enabled.



- Plot Statistics:** By default, all gates show the **% Gated** statistic. The selection persists on a per Workspace basis.  
 When a statistic option is selected, all plots with gates are updated to show the selected statistic. Selecting a new statistic unselects the previously selected one.
- Statistics tables:** Individual **Plot statistics** and **Workspace statistics** can be selected at the same time. When multiple statistics tables are selected, the selection and deselection of statistics in the **Statistics** tab applies to all tables.  
 Where a statistic is selected in one statistics table and not in another, the checkbox is in the indeterminate state.  
 The selection or deselection of the **Experiment name**, **Group name**, **Sample name**, and **Plot title** also updates the selections in the **Customize statistics** panel.

### Behavior in Results view

- When the **Results** view (Chapter 7, “Sample List view”) is in focus, all statistics can be selected and unselected. The **Results** group is only available when the Results view is in focus.



- A statistical value is displayed in the **Results table** when the associated checkbox is checked in the **Statistics** tab. Unselecting the checkbox removes the entire column from the **Results table**.

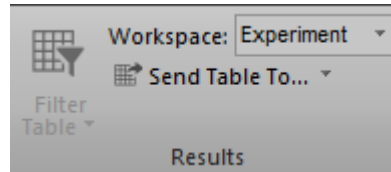
### Persistence of selections

The persistence of selections in the Statistics tab is context dependent.

- For Workspace selections, the selections for on-plot statistics and for each statistics box persist on a Workspace level.
- For the Results view, all selections (including the filter selection) persist on an Experiment-level.

## Results group

The **Results** group contains the controls to turn the **Results table filter ON** and **OFF**, to edit the filter criteria, to select the Workspace on which the **Results table** is based, and to select where to send the **Results table** (“Results table” on page 229).

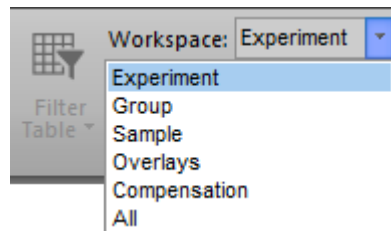


The **Results** group is only visible when the **Results** view is in focus.

### Workspace filter dropdown

The **Workspace filter** dropdown enables you to select the **Workspace** on which the **Results table** is based.

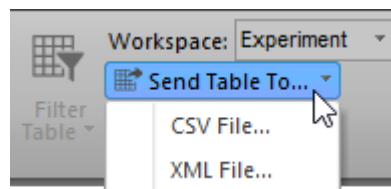
The options include **Experiment**, **Group**, **Sample**, **Overlays**, **Compensation**, and **All**. When **All** is selected, all options are displayed. By default, **Experiment** is selected.



### Send Table To...dropdown

The **Send Table To...** function exports all results that are displayed in the **Results** view, including all relevant statistics for the region and gate type, and filters; grouping is not included.

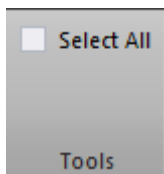
The **Send Table To...** dropdown enables you to export the results from the experiment as a **CSV File** or as an **XML File**.



By default, all experimental metadata, statistics, and keywords are exported as an XML file.

## Tools group

The **Tools** group lets you enable all statistics with a single click.



**Select All:** Enables all statistics. The **Select All** option is enabled only when the **Results** view is in focus or if a Statistics box is selected on the **Reports** view or the **Workspace** view. The checkbox is disabled for all other instances.

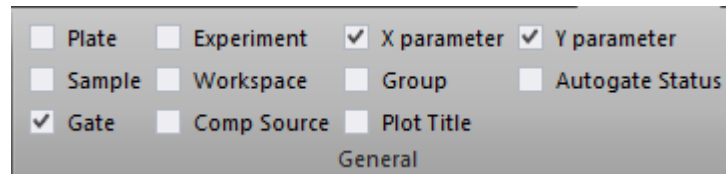
When **Select All** is selected, then all statistics are selected. If any statistics are deselected, the **Select All** option becomes deselected.

Deselecting the **Select All** option deselects all statistics.



## General group

The statistics in the **General** group, when selected, are displayed on plots and in the **Results table**.

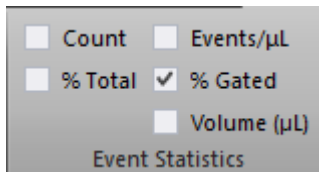


If a statistics table is not selected, this group is disabled on the Workspace view.

- **Plate:** Displays the Plate name (enabled for Plate Experiments only)
- **Experiment:** Displays the Experiment name.
- **Group:** Displays the Group name.
- **Sample:** Displays the Sample name.
- **Workspace:** Displays the Workspace from which the statistics is calculated.
- **Comp Source:** Displays the Compensation Source for the Results, either calculated in the Experiment or embedded with FCS files
- **Plot Title:** Displays the plot title for all gates.
- **Gate:** Displays the name of the gate for which the statistics is calculated.
- **X Parameter:** Displays the X parameter name as shown on the plot that contains the gate.
- **Y Parameter:** Displays the Y parameter name as shown on the plot that contains the gate. If the Y parameter is from a Histogram plot, **Count** or **% of max** is displayed as the parameter name, depending on what is selected for that plot.
- **Autogate Status:** Displays the status of an autogate.

## Event Statistics group

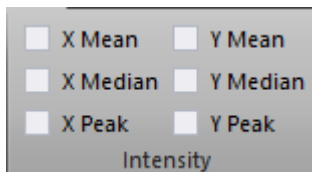
The statistics in the **Event Statistics** group, when selected, are displayed on plots and in the **Results table**.



- **Count:** Displays the number of events in gates.
- **% Total:** Displays the percentage of total events.
- **Events/μL (Concentration):** Displays the number of events per μL.
- **% Gated:** Displays the percentage of events within the gate.
- **Volume (μL):** Displays the volume of sample from which the data are derived.

## Intensity group

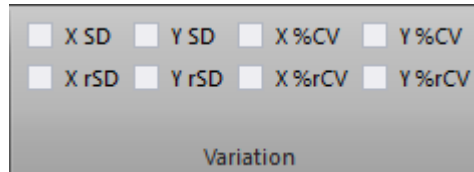
The statistics in the **Intensity** group, when selected, are displayed on plots and in the **Results table**.



- **X Mean:** Displays the arithmetic mean for the X-axis of a gate.
- **Y Mean:** Displays the arithmetic mean for the Y-axis of a gate. When the gate is on a single-parameter plot, the entry on the **Results table** is **N/A**.
- **X Median:** Displays the median for the X-axis of the gate.
- **Y Median:** Displays the median for the Y-axis of the gate. When the gate is on a single-parameter plot the entry on the **Results table** is **N/A**.
- **X Peak:** Displays the peak channel (mode) for the X-axis of the gate.
- **Y Peak:** Displays the peak channel (mode) for the Y-axis of the gate. When the gate is on a single-parameter plot, the entry on the **Results table** is **N/A**.

## Variation group

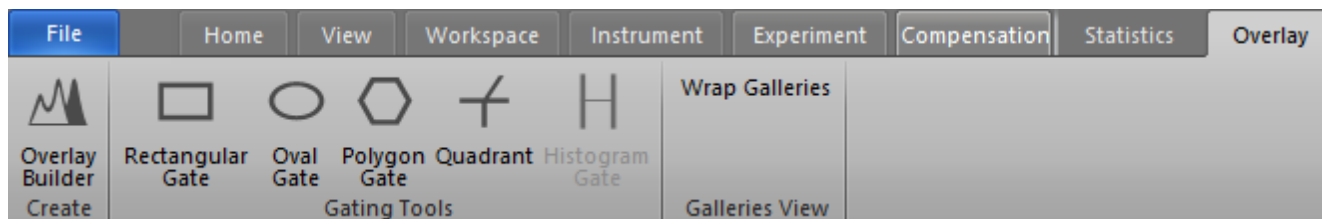
The statistics in the **Variation** group, when selected, are displayed on plots and in the **Results table**.



- **X SD**: Displays the standard deviation for the X-axis of a gate.
- **Y SD**: Displays the standard deviation for the Y-axis of a gate. When the gate is on a single-parameter plot, the entry on the **Results table** is **N/A**.
- **X %CV**: Displays the %CV (percentage coefficient of variation) statistics for the X-axis of a gate.
- **Y %CV**: Displays the %CV (percentage coefficient of variation) statistics for the Y-axis of a gate. When the gate is on a single-parameter plot, the entry on the **Results table** is **N/A**.
- **X rSD**: Displays the robust standard deviation statistics for the X-axis of a gate.
- **Y rSD**: Displays the robust standard deviation statistics for the Y-axis of a gate. When the gate is on a single-parameter plot, the entry on the **Results table** is **N/A**.
- **X %rCV**: Displays the percentage robust CV (coefficient of variation) statistics for the X-axis of a gate.
- **Y %rCV**: Displays the percentage robust CV (coefficient of variation) statistics for the Y-axis of a gate. When the gate is on a single-parameter plot, the entry on the **Results table** is **N/A**.

## Overlay tab

**Overlay** ribbon tab contains the tools used for creating **Overlays**, which are used for comparing data by superimposing selected plots. Each **Overlay** consists of an **Overlay plot** and an associated **Overlay gallery** (Chapter 9, “Overlays”).



**Overlay** tab contains three functional groups:

- **Create** (“Create group” on page 74)
- **Gating Tools** (“Gating Tools group” on page 111)
- **Galleries View** (“Galleries View group” on page 111)

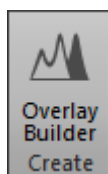
---

**Note:** The **Overlay plots** and the associated **Overlay galleries** are displayed in the **Overlay view** (Chapter 9, “Overlays”).

---

### Create group

**Create** group contains the **Overlay Builder** button.



- **Overlay Builder** launches the **Overlay Builder dialog** (“Overlay Builder” on page 260), which lets you select the **Samples** and **Plots** for inclusion in an **Overlay plot**.
- **Overlay Builder** is only enabled when at least one plot exists on the Experiment Workspace and the instrument is not acquiring. It is disabled when a Compensation control is active.

## Gating Tools group

**Gating Tools** group contains the controls used for inserting **Rectangle**, **Oval**, **Polygon**, **Quadrant**, and **Histogram gates** into **Overlay plots**. These controls are only enabled if at least one **Overlay plot** exists.



- Rectangular Gate:** Enables you to draw a **Rectangular gate** on a dual-parameter plot.
- Oval Gate:** Enables you to draw an **Oval gate** on a dual-parameter plot.
- Polygon Gate:** Enables you to draw a **Polygon gate** on a dual-parameter plot.
- Quadrant Gate:** Enables you to draw a **Quadrant gate** on a dual-parameter plot.
- Histogram Gate:** Enables you to draw a **Histogram gate** on a Histogram plot.

---

**Note:** The controls in the **Gating Tools** group are highlighted when selected. They revert to their standard color when you click the **Overlay plot** to create a gate, select another control, or press the **ESC** key.

---

## Galleries View group

**Galleries View** group contains the **Wrap Galleries** button.

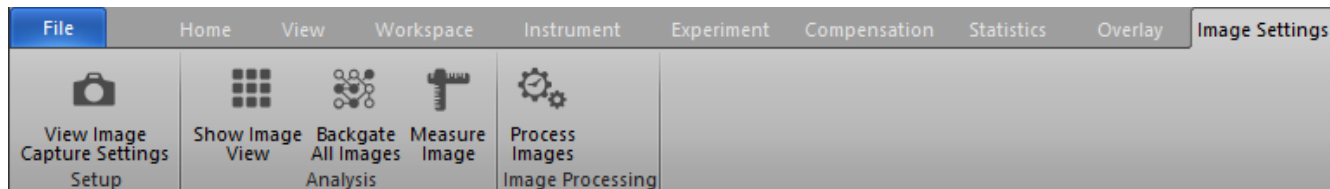


When **Wrap Galleries** is selected, the plots in the **Overlays view** all fit within the width of the **Galleries** area. If there are too many plots to be shown vertically, vertical scroll bars are shown.

When the gallery is wrapped, the **Wrap Galleries** button is highlighted in blue. The button is only enabled if the **Gallery plot view** or the **Overlay view** is in focus.

## Image Settings tab

**Image Settings** ribbon tab provides options to set up cell image capture and acquisition settings, view captured images, enable image backgating, measure image dimensions, and process images.



The **Image Settings** tab is only visible when the Attune™ desktop is in view and one of the following criteria are met:

- An Attune™ CytPix™ Flow Cytometer is in use or is selected as the active instrument type in **Configuration** options (“Hardware/Virtual laser configuration” on page 696)
- When an Experiment that was created with an Attune™ CytPix™ Flow Cytometer is active.

The **Image Settings** tab is not visible on the **Main Menu**, **Login screen**, **Performance Test views**, or **Filter Configuration views**.

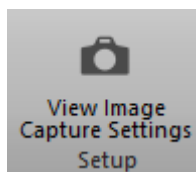
The **Image Settings** tab is not available for the Attune™ NxT Flow Cytometer (unless viewing an Experiment that was created using an Attune™ CytPix™ Flow Cytometer).

The **Image Settings** tab contains three functional groups:

- **Setup** (“Setup group” on page 112)
- **Analysis** (“Analysis group” on page 113)
- **Image Processing** (“Image Processing group” on page 113)

### Setup group

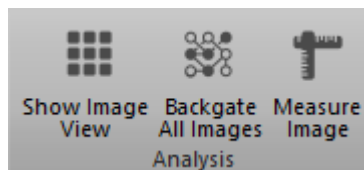
The **Setup** group contains the **View Image Capture Settings** button.



- **View Image Capture Settings** toggles the display of the **Image Capture Settings** panel (Chapter 16, “Image Capture Settings panel”), which lets you adjust settings that affect the capture of images.
- When the **Image Capture Settings** panel is displayed, the **View Image Capture Settings** button is in the **ON** state (blue).  
When the panel is closed, the **View Image Capture Settings** button is in the **OFF** state.
- By default, when the instrument model is set to Attune™ CytPix™ Flow Cytometer and the first Experiment is activated, the **View Image Capture Settings** button is in the **ON** state. After this, the **ON/OFF** state of the button persists with user settings.

## Analysis group

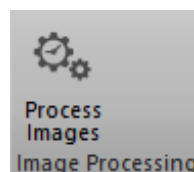
The **Analysis** group lets you show and hide the **Image View**, to backgate images, and to measure images.



- **Show Image View** toggles the display of the **Image View** tab (Chapter 10, “Image View”).  
When the **Image View** tab is displayed, the **Show Image View** button is in the **ON** state (blue).  
When the **Image View** tab is closed, the **Show Image View** button is in the **OFF** state.  
By default, when the instrument model is set to Attune™ CytPix™ Flow Cytometer and the first Experiment is activated, the **Show Image View** button is in the **ON** state. After this, the **ON/OFF** state of the button persists with user settings.
- **Backgate All Images** toggles the backgating of all imaged events on Workspace plots. Backgating enables you to identify the events on the plot in Workspace view that are associated with the captured cell images.  
By default, backgated events appear on the associated plot in red. For more information, see “Image backgating” (“Image backgating” on page 179).  
When image backgating is applied, the **Backgate All Images** button is in the **ON** state (blue). Otherwise, the button is displayed in its **OFF** state.  
Image backgating persists at the Workspace level for an Experiment, Group, or Sample.  
The **Backgate All Images** button is enabled only if the selected Sample’s listmode specification contains the **ImageFlag** and **Event** parameters, and the Workspace is in focus/active.
- **Measure Image** enables you to insert a measurement ellipse on the active image in the **Image view** to measure the area of the cell in the image (“Measure image tool” on page 280).  
When using the measurement tool to insert the measuring ellipse, the button is displayed in the **ON** state (blue).

## Image Processing group

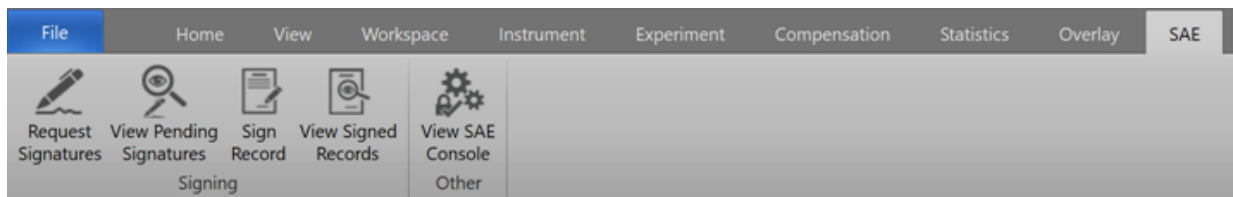
**Image Processing** group contains the **Process Images** button, which opens the **Process Images** dialog (Chapter 23, “Process Images dialog”).



The **Process Images** dialog is used to process captured images from selected **Samples** and **Populations** using system-provided **Image Processing Models**.

## SAE tab

The **SAE** ribbon tab provides tools to manage signatures (e-Signing) and to access the **SAE Administrator Console** (“SAE Administrator Console” on page 880), where you can configure the Security, Auditing, and e-Signature (SAE) settings. **SAE** tab is available only when an SAE user is signed in.



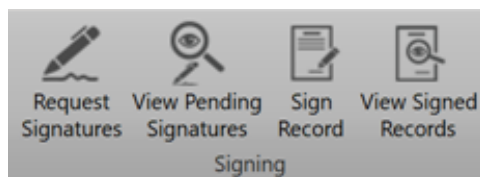
The **SAE** tab is divided into two functional groups:

- **Signing** group (“Signing group” on page 114)
- **Other** group (“Other group” on page 115)

The **SAE** tab is only visible on the **Main Menu** and **Workspace views**. During acquisition, when the plate is paused, or when in automation mode, all buttons are disabled.

### Signing group

The **Signing** group enables you to request e-Signatures for the active Experiment, view pending signature requests, sign the active Experiment, and view signed records in the **e-Signature Record Report History**.



**Request Signatures:** Opens the **Request Signatures** dialog (“Request Signature dialog” on page 806), which enables you to request signatures for the active Experiment. This button is only enabled if **e-Signatures** are enabled from the **SAE Administrator Console**, actions are set up that require signatures, and an Experiment is active.

**View Pending Signatures:** Opens the **Sign Records** dialog (“Sign Records dialog” on page 808), which enables you to view and sign pending signature requests. This button is only enabled if **e-Signatures** are enabled from the **SAE Administrator Console**.

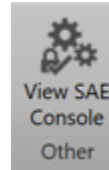
**Sign Record:** Opens the **Sign Experiment** dialog (“Sign Experiment dialog” on page 810), which lets you sign the active Experiment. This button is only enabled if **e-Signatures** are enabled from the **SAE Administrator Console** and an Experiment is active.

**View Signed Record:** Opens the **e-Signature Record Report History** dialog (“e-Signature Record Report History dialog” on page 812), which lets you view and print signed records.



## Other group

The **Other** group contains the **View SAE Console** button, which opens the **SAE Administrator Console** to enable you to configure **Security, Auditing, and e-Signature settings**, and to view auditing and signature records.

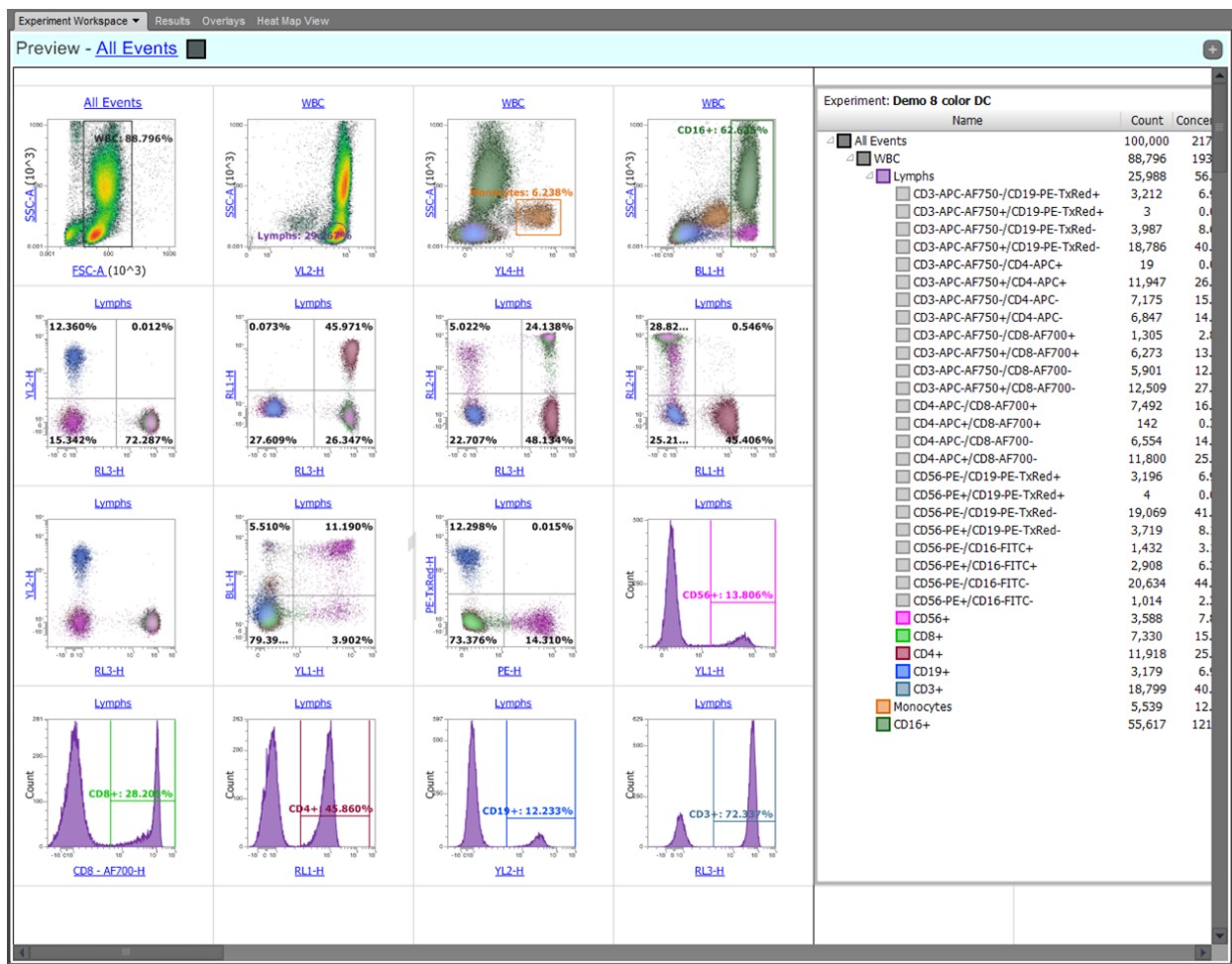


**View SAE Console:** Opens the **SAE Administrator Console** (“SAE Administrator Console” on page 880) using the system’s default browser.

The **SAE Administrator Console** enables those with SAE administrative role to configure **Security, Auditing, and e-Signature settings**, to view application action records, application object records, and signature history, and to set up advanced settings including email notification, LDAP user management synchronization, and SAE archive settings.

## Overview

The **Workspace view** displays the plots, gates, and statistics associated with the current Sample, as well as text fields and image files. It also contains the **Preview panel** (“Previews” on page 132), which is displayed by default as a minimized bar.



The Workspace view contains these elements:

- **Workspace** (“Workspace selection” on page 118)
- **Workspace objects** (“Workspace objects” on page 120)
- **Previews** (“Previews” on page 132)
- **Plots** (“Plots” on page 138)

- **Gates** (“Gates” on page 145)
- **Workspace statistics** (“Workspace statistics” on page 175)
- **Workspace images and text** (“Workspace images and text” on page 178)
- **Cell image container** (“Cell image container” on page 181)

You can access **Workspace view** and **Workspace object context menus** when you right-click any empty area (i.e., white space) in the **Workspace view** or when you right-click **Workspace objects**.

## Workspace behavior

- The **Workspace view** is always the first tab of the **Main Application area** (“Main application workspace” on page 56). It cannot be undocked from the main application and the associated ribbon bar (see “Docking locations” on page 58).
- The **Workspace view** can show one of three independent workspaces, the **Sample Workspace**, the **Group Workspace**, or the **Experiment Workspace** (“Workspace levels” on page 118)
- The **Workspace** can be displayed in two different modes, **Freeform** or **Auto Layout** (“Auto Layout mode” on page 127). These enable automatic or user-specified placement of objects on the **Workspace** depending on the mode selected.
- The **Workspace** acts as a single worksheet that is split into “pages” when printed. The page breaks can be visualized by selecting **Print Area** in the **View tab** (“View tab” on page 77).
- Changes to the **Workspace** are saved automatically if the active **Workspace** changes, if the program is exited, if the active Sample changes, before printing, or if one minute has elapsed since the last Workspace save action.
- When a Compensation sample is active, all Workspace functionality is disabled except as specified in Chapter 21, “Compensation”.
- The **Workspace view** always moves to the front of any docked panels at the beginning of acquisition for each sample. The **Experiment Workspace** becomes active and the dropdown menu items become disabled.

---

**Note:** The **Preview panel** shows previews of all possible permutations of the Histogram and Precedence Density plots based on the parameters selected in the **Workspace** tab, which enable you to quickly add these plots to the **Workspace** as described in “Previews” on page 132.

---

## Workspace selection

### Workspace levels

The **Workspace view** can show one of three independent workspaces, the **Sample Workspace**, the **Group Workspace**, or the **Experiment Workspace**.

- **Experiment Workspace** is a Workspace that is common to all members of the current Experiment.
- **Group Workspace**, when created, is common to all members of the current Group.
- **Sample Workspace** is unique to each Sample.

### Workspace view label

The **Workspace view** tab on the **Main Application** area is labeled according to the level of the **Workspace** displayed: **Sample Workspace**, **Group Workspace**, or **Experiment Workspace**.

- When a Compensation sample is open, the Workspace is labeled **Compensation Workspace**, and the dropdown selection menu is not present.

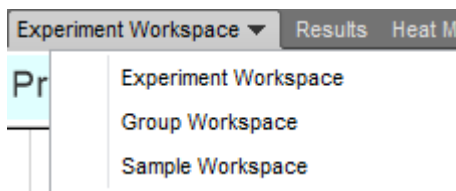
Compensation Workspace

- By default, the **Experiment Workspace** is shown when the Sample is first opened. **Group Workspaces** and **Sample Workspaces** are created upon first use.

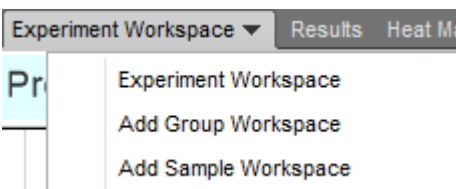
Experiment Workspace ▼

### Workspace selection menu

- The **Workspace selection dropdown** on the **Workspace view** tab lets you select the level of **Workspace** displayed:



- **Experiment Workspace**
- **Group Workspace**
- **Sample Workspace**
- If the **Group** or **Sample Workspaces** do not exist, the workspace label is displayed as **Add Group Workspace** or **Add Sample Workspace**.



- If you select a **Group Workspace**, that **Workspace** is displayed. If a **Group Workspace** does not exist for the current Group, a default **Group-level Workspace** is created.
- If you select the **Sample Workspace** option, the **Sample Workspace** for that Sample is displayed. If a **Sample Workspace** does not exist for the current Sample, a default **Sample-level Workspace** is created.
- If you select the **Experiment Workspace**, the current **Experiment Workspace** is displayed.
- When a new **Group Workspace** is created, the **Experiment Workspace** is used as a template, and all plots, gates, and statistics are reproduced.
- When a new **Sample Workspace** is created, the **Group Workspace** is used as a template. If the **Group Workspace** does not exist, then the **Sample Workspace** uses the **Experiment Workspace** as a template.
- When a new **Group** or **Sample Workspace** is created, an indicator for the Group or Sample Workspace is displayed in the **Experiment Explorer** as described in Chapter 11, “Experiment Explorer”.

## Workspace objects

### Create Workspace objects

- The **Workspace** can contain plots, gates, statistics, text boxes, and images (i.e., **Workspace objects**), all of which can be created using the options available on the **Workspace ribbon tab** (“Workspace tab” on page 81).

If you are using an Attune™ CytPix™ Flow Cytometer or the current Experiment supports imaging, you can add **Cell Image Containers** to the Workspace, which act like plots and let you add cell images specific to the Sample (“Cell image container” on page 181).

You can use the **Workspace view context menu** (“Workspace view context menu” on page 185) to add the Workspace objects onto the Workspace.

- The **Plot context menu** (“Plot context menu” on page 189) enables you to add plot-specific statistics.
- The **Preview panel** enables you to add plots to the Workspace (“Add Plots to Workspace – Single Plot selection” on page 134).
- In the **Freeform** mode (“Freeform mode” on page 126), you can add Workspace objects (i.e., plots, text boxes, images, or statistics) of any size to the Workspace. These objects can be freely resized and moved on the Workspace. There is no automatic arrangement or alignment of objects.
- In the **Auto Layout** mode (“Auto Layout mode” on page 127), the Workspace is split into grid slots. By default, new Workspace objects are placed into the next available grid space and are sized to one grid square.
- Single plots that are added using the **Previews** option (“Add Plots to Workspace – Single Plot selection” on page 134) or the **daughter plot** option (“Create daughter plot” on page 195) are inserted at the next available space of sufficient size after the last workspace object. Objects are created using the default object size.

## Select Workspace objects

- To select a **Workspace objects**, click within the **gray bounding rectangle** that is displayed when you move the mouse pointer over the object.

When you select a **Workspace object**, it is displayed with a **light blue selection rectangle**.

Workspace objects that are not selected do not have a selection rectangle displayed.

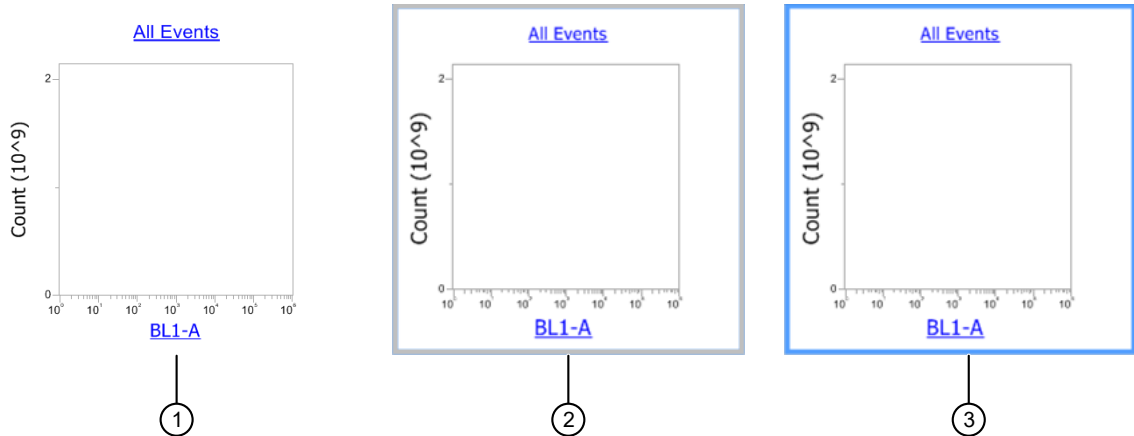
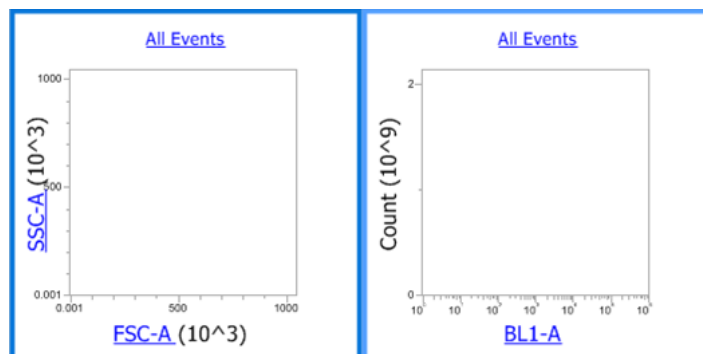
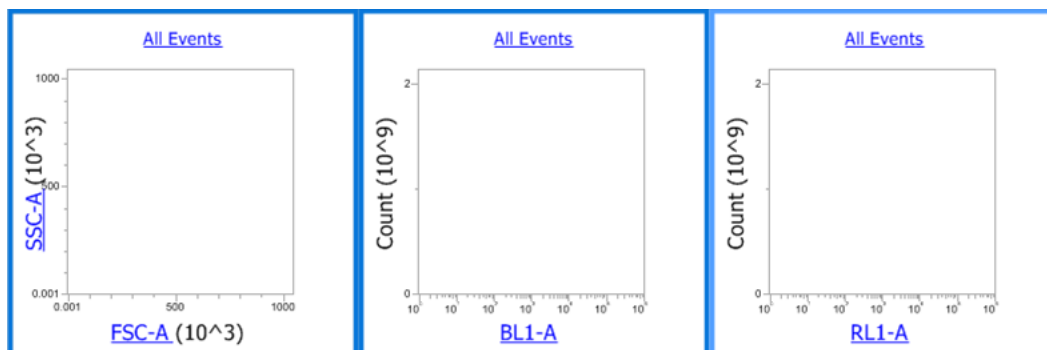


Figure 13 Selecting Workspace objects

- ① Workspace object (Plot) that is not selected.
  - ② Workspace object shows the gray bounding rectangle when moused over.
  - ③ Workspace object shows the light blue selection rectangle when selected.
- To select multiple objects in the Workspace, click the desired **Workspace objects** while holding down the **Ctrl** key. Alternatively, you can drag a **selection rectangle** around several objects with the mouse. **Ctrl+A** key combination on the keyboard selects all Workspace objects on the current Workspace. Objects which are part of the multiselect are bounded by **dark blue multiselect rectangles**.



- When multiple objects are selected (either individually clicked or selected by dragging the selection rectangle), the last selected object remains light blue to designate it as the **reference item for align options**.



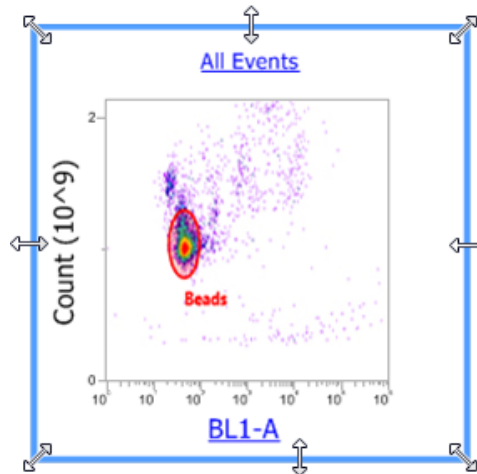


## Resize Workspace objects

- To resize single or multiple **Workspace objects**, select the object, then move the mouse pointer over the edge of the bounding rectangle.

**Vertical**, **horizontal**, and **diagonal resize arrows** are displayed as you move the mouse pointer over the appropriate area.

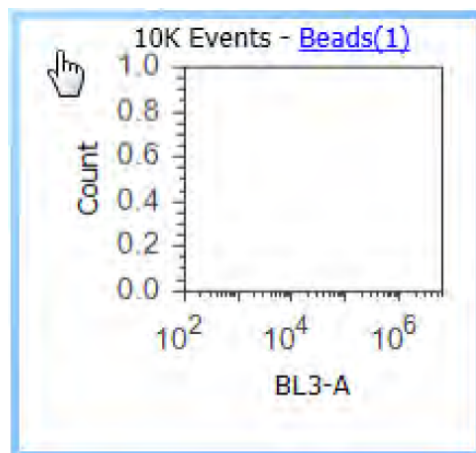
The **resize arrows** are also displayed when the mouse pointer is hovered over the bounding rectangle of non-selected objects; when clicked, this selects the object and begins resizing.



- When the resize arrows are displayed, the object can be dragged to resize in the direction indicated by the arrow. All selected objects are resized in the same dimension and magnitude.
- Dual parameter plots and images always maintain their aspect ratio when resized. Histograms can be stretched horizontally. They can be stretched vertically, but the width of the object must always be greater than or equal to its height. Text fields and statistics can be stretched in any direction and any ratio.
- Pressing the **Shift** key and dragging the diagonal resizing arrows resizes the selected object and maintains its aspect ratio.
- When multiple Workspace objects are selected, resizing options are determined by the object being manipulated. All other objects in the selection are bound by the resizing constraints described above.
- If mixed object types are selected when manipulating an object whose aspect ratio is not limited, histogram and dual parameter plots are resized with the aspect ratio constraints described above. Where necessary, requirements for width take precedence over height to ensure that the resizing constraints are met.

## Move Workspace objects

- Moving the pointer over a **Workspace object** displays the **select cursor** in the shape of a pointing finger, except when the pointer is over a line or the select point of a gate. In such cases, **resize**, **move**, or **rotate cursors** are displayed as described on “Move and resize gates” on page 148.
- When the **select cursor** is displayed, you can click and drag single or multiple selected **Workspace objects** to a new position.



## Delete, format, or duplicate Workspace objects

### Delete Workspace objects

- To delete selected **Workspace objects**, press the **Delete** key on the keyboard.
- Alternatively, right-click the object to open the **Object context menu**, then select the **Delete** option.

### Format Workspace objects

- You can format the selected **Workspace object** using the **Customize panel** options (Chapter 15, “Customize panel”).
- Alternatively, right-click the object to open the **Object context menu**, then select the desired formatting option.

### Duplicate Workspace objects

- To duplicate **Workspace objects**, select one or more objects, then drag them as you press and hold the **Ctrl** key on the keyboard.
- Objects can also be duplicated using the keyboard combinations **Ctrl+C** (copy) followed by **Ctrl+V** (paste).
- Alternatively, copy the object using the **Copy** command on the object’s **Object context menu**, then paste it using the keyboard combination **Ctrl+V**.

## Paste Workspace objects

- Use the **Ctrl+V** keyboard combination to paste the object at same location as the original. As each new object is pasted, it is inserted cascading downwards and to the right of the original location.
- Objects can be cut or copied from any Workspace, then pasted to any other Workspace.
- Plots are pasted with any gates present. If a gated plot is copied to a new location where the parent gate does not exist, the pasted plot is not be gated and displays all events.
- If a Plot statistics box is pasted to a new Workspace and the target Workspace does not contain a copy of the original plot, the statistics box updates to show all gates present in the current Workspace. The Plot axis and Gate fields are removed from the statistics box header (if originally present).
- Pasted objects are selected by default after the paste action has completed.

## Workspace view layout

### Resize the Workspace view

There are two ways to resize the **Workspace view**:

- When the mouse cursor is over the **Workspace** or when the Workspace is in focus, turn the **mouse scroll wheel** as you hold down the **Ctrl** key to zoom in or out.
- Use the **Size slider** on the **Application status bar** (“Size slider” on page 70) to resize the **Workspace view** when the **Workspace** is in focus.

### Workspace view modes

The **Workspace** can be displayed in two different modes, **Freeform** or **Auto Layout**. These modes enable automatic or user-specified placement of objects on the **Workspace** depending on the mode selected.

The **Workspace view mode** is controlled by the **Layout** buttons available on the **Workspace ribbon tab** (“Grid options” on page 87).

### Freeform mode

**Freeform** mode lets you add plots, text boxes, images, or statistics of any size anywhere on the **Workspace**. These objects can be freely resized and moved on the **Workspace**. There is no automatic arrangement or alignment of objects.

- To add a custom-size object to the **Workspace**, select the desired object type from the **Workspace ribbon tab** (“View tab” on page 77) and then draw the area in the **Workspace** where you want to insert the selected object. When you let go of the mouse button, the object is inserted into the area drawn.
- Selecting an object type and then single-clicking on the **Workspace** places the default-sized object at the position clicked.

The default size for new plots, statistics boxes, and text boxes in the **Freeform** mode is 240 × 240 pixels with objects arranged in 3 columns × 4 rows per page.

- In the **Freeform** mode, objects can be moved and positioned anywhere in the **Workspace**, and they can be positioned across multiple pages.
- You cannot print a **Workspace** in the **Freeform** mode.

## Auto Layout mode

In the **Auto Layout** mode, the **Workspace** is split into grid slots, which contain the **Workspace objects** of fixed size and placement. **Auto Layout** is the default option.

### Layout behavior

- Only one object can appear in each grid slot. By default, new **Workspace objects** are placed into the next available grid space and are sized to one grid square.
- When switching from **Freeform** to **Auto Layout** mode, any objects present on the **Workspace** are placed into grid positions in the following order:
  - Plots based in Workspace insertion order
  - Plots statistics based on Plot insertion order
  - Workspace statistics
  - Text boxes
  - Images
- When switching from **Auto Layout** to **Freeform** mode, objects remain in their current position.
- The **Workspace** is initially setup as one page view. More pages are added automatically as existing pages are filled and as the **Workspace zoom** is changed.

### Grid size

- The number of available grid slots per Workspace page is defined by the **Grid Size** options on the **Workspace ribbon tab** (“Grid options” on page 87).
- By default, the grid size is set to 12 grid slots per page.
- You can select a larger grid size allowing only 4 items per page, or a smaller grid size allowing up to 20 items per page at the default sizes.

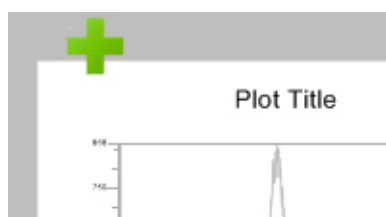
### Object size

- By default, objects on the **Workspace** are sized to one grid square when the **Workspace view** is in the **Auto Layout** mode.
- Objects on the **Workspace** can be resized, but they have limited resize options. When an object is resized, the top left corner of the object is pinned to its current position. Objects cannot be resized across pages or resized to occupy more than one page.

- The resize options available for various **Workspace objects** are listed below, where the sizes are defined as (Rows × Columns).
  - **Dual parameter plots** are constrained to remain square and can be resized to take occupy 1, 4 (2 × 2), 9 (3 × 3) or 16 (4 × 4) squares.
  - **Histogram plots** are either constrained to remain square and occupy 1, 4 (2 × 2), 9 (3 × 3), or 16 (4 × 4) grid spaces, or they can be made rectangular and occupy 2 (1 × 2), 3 (1 × 3), 4 (1 × 4) or 2 × 2), 6 (2 × 3), 8 (2 × 4), 12 (3 × 4), or 16 (4 × 4) squares.
  - **Statistics boxes** and **images** can be resized up to the maximum number of grids (5 × 4) per page. For images, the aspect ratio is maintained in the grid containers.
- When increasing grid sizes (from more slots per page to less slots per page) or changing page orientation, any object that exceeds the allotted slots in the horizontal and/or vertical directions is resized to 1 × 1.

### Moving objects in Auto Layout mode

- Object arrangement precedence is dictated by grid position and page number. The position with the highest precedence is the top left corner of the first page, and the precedence decreases from left to right, top to bottom.  
The grid position of objects larger than 1 × 1 is dictated by the top, leftmost position occupied by that object.

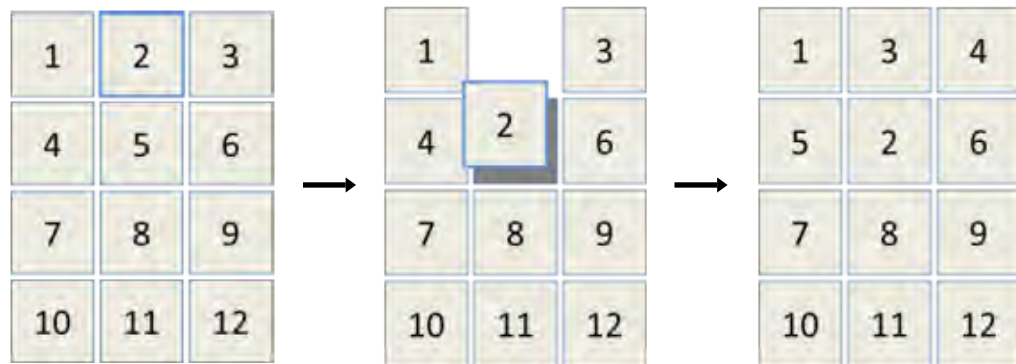


- When moving an object, a **gray box** indicates the grid location into which the object will be inserted or expanded.
- When multiple objects are selected, only the object with the highest precedence is displayed, and a **plus icon** indicates that multiple objects have been selected.
- If the objects have been selected using the **Ctrl+click** method, then the first selected object is displayed when dragging them.  
If the objects have been selected by dragging a selection area, then the object with the highest precedence based on Workspace location (top to bottom, left to right) is displayed.
- After objects have been resized or moved, the selection indicators are maintained on the selected objects.

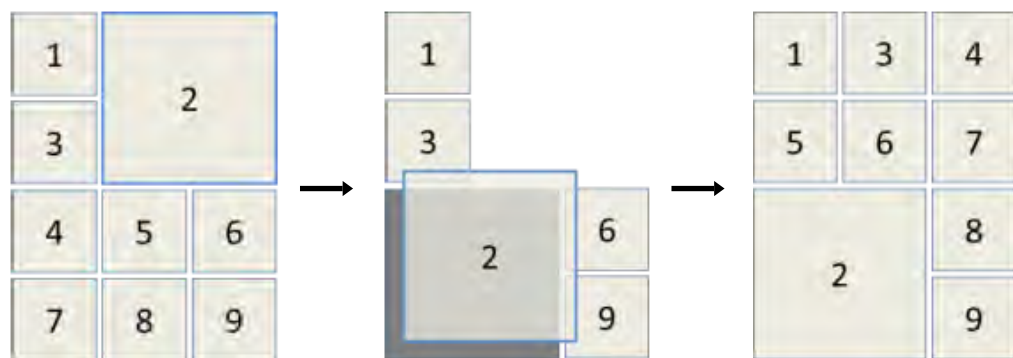
## Rules for rearranging objects in Auto Layout mode

When a resized object causes another object to be displaced or a deleted object results in a blank grid slot, the objects on the **Workspace** are rearranged according to the rules listed below.

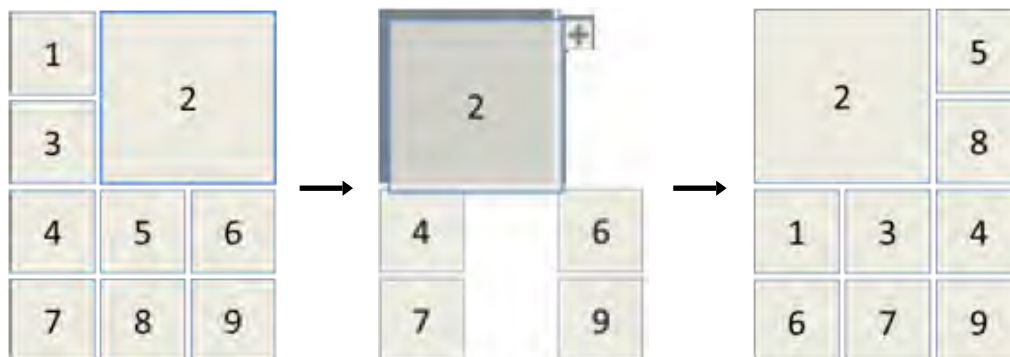
1. The displacing objects are always inserted in the position indicated by the gray box.
  2. Displaced objects are always placed in the highest precedence available grid position.
  3. The size of the object cannot span more than one page
  4. Objects will always fill blank positions whenever possible.
- **Example 1:** Plot 2 is a dot plot and it is dragged over Plot 5. Plot 3 moves to fill position 2, Plot 4 moves to fill position 3, and Plot 5 moves to fill position 4. Plot 2 takes the position of Plot 5.



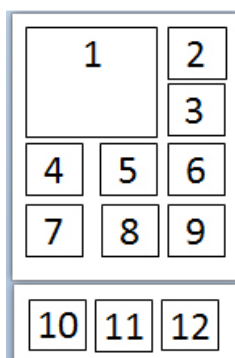
- **Example 2:** Plot 2 is a dot plot occupying 4 grid positions and it is dragged over Plot 4. Plots 3, 4, 5, 6, 7, 8, and 9 move to refill the displaced space.



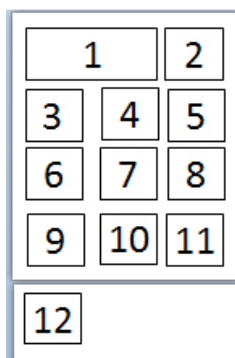
- Example 3:** Plot 2 (a dot plot occupying 4 grid positions) along with Plots 5 and 8 are dragged over position 1. Plots 1, 3, 4, 6, 7, and 9 move to refill the displaced spaces.



- Example 4:** Plot 1 is a dot plot and it is resized to be 4 grid slots. Plots 2 and 3 automatically rearrange to optimally fill the space. Plots 10–12 are pushed onto the next page, which appears below the current page.

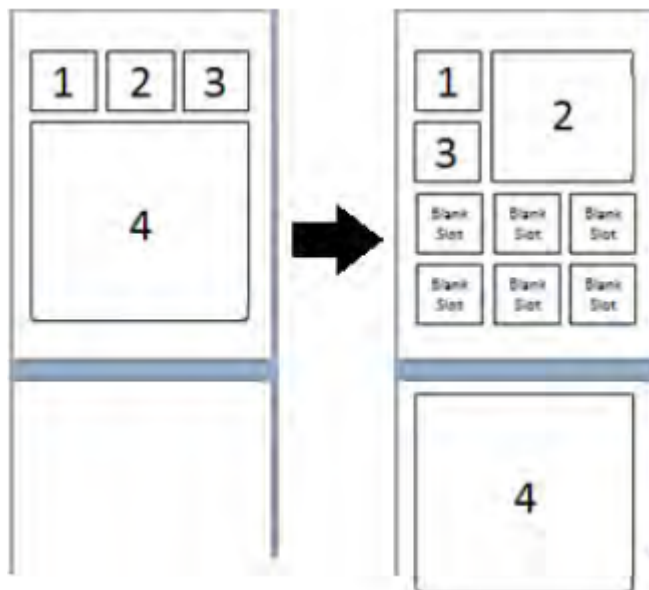


- Example 5:** Plot 1 is a histogram and it is resized to 2 grid slots. Plots automatically rearrange as required, and Plot 12 is pushed onto the next page, which appears below the current page.





- **Example 6:** Plot 2 is dot plot and it is resized to be 4 grid slots. Plots automatically rearrange as required, and Plot 4 is pushed onto the next page, which appears below the current page. Because there are no other plots available to occupy them, blank slots are left on the original page



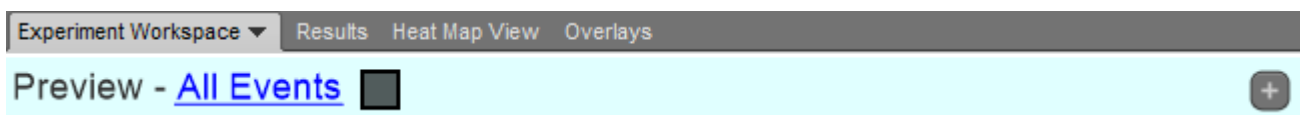
## Previews

### Overview

The **Preview panel** of the **Workspace** displays all permutations of Histogram (“Histogram plot” on page 139) and Precedence Density plots (“Precedence density plot” on page 142) based on the parameters selected in the **Workspace ribbon tab** (“Workspace tab” on page 81). This provides an easy way of adding plots to the current **Workspace**.

### Access the Preview panel

The **Preview panel** is at the top portion of the **Workspace view**. By default, the panel is displayed as a minimized bar.



- When opened, the **Preview panel** occupies 50% of the vertical space of the **Workspace view**. The size of the panel cannot be changed.
- The **Preview panel** cannot be opened during acquisition. The panel closes when **Tube acquisition** or **Plate acquisition** buttons are selected on the **Collection panel** (Chapter 12, “Collection panel”).
- The **Preview panel** is disabled in the **Compensation Workspace**.
- When the **Preview panel** is minimized, the **Expand** button is shown in the collapsed panel.



The **Preview panel** can be expanded by clicking anywhere, except for the **gate hyperlink** on the **title bar** of the collapsed **Preview panel**. The expanded setting of the **Preview panel** does not persist between user sessions.

- When the **Preview panel** is expanded, the **Collapse** button is shown on the panel.




The **Preview panel** be collapsed by clicking anywhere on the **title bar**, except for the **gate hyperlink** of the expanded **Preview panel** or when the panel is closed.

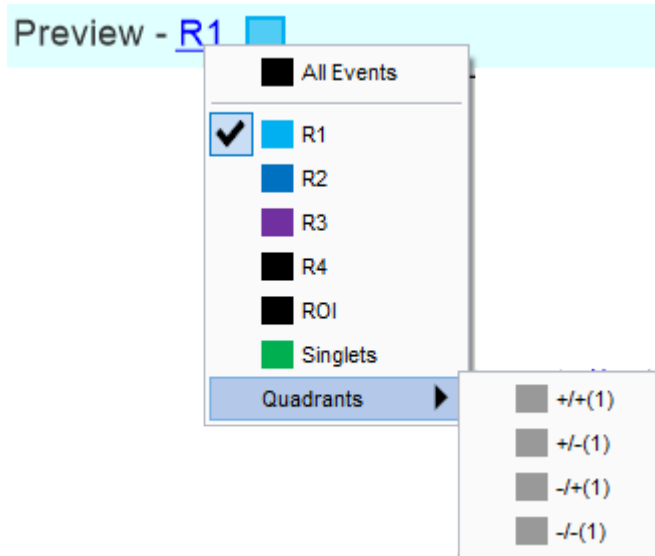
- When the **Preview panel** is disabled during acquisition or in the **Compensation Workspace**, the **gate hyperlink** and the **Expand/Collapse** buttons are also disabled.

## Preview panel title bar

- The **Preview panel title bar** displays “Previews – <Gate Name> <Gate Color>” where <Gate Name> is the name of the selected gate used to filter the **Preview plots**. The **Gate Name** is a hyperlink control.

Preview - [Singlets](#) 

- Clicking on the **Gate Name** or the **Gate Color** on the **Preview panel title bar** opens a menu of all available gates and an option for **All Events** at the top of the list. Quadrant gates are shown in a submenu under the title **Quadrants**.



## Preview plots

### Layout of Preview plots

- **Preview plots** are arranged in a grid organized such that all plots share the same X parameter within a column and the same Y parameter within a row.



- The single-parameter **Histogram plots** are displayed in the first row.
- **Dual-parameter plots** show all possible combinations of available parameters.
- The plot order is based on parameter order, and starts with the first parameter in Column 1, Row 1 and ends with the Nth parameter in Column N, Row N.
- **Self vs. Self dual parameter plots** are not shown in the **Preview panel**. They are represented by white spaces that form a diagonal from the top left to the lower right of the panel.
- The default scale for **Preview plots** is set to **logarithmic**.

## Resize Preview plots

Turning the **mouse scroll wheel** when holding down the **Ctrl** key resizes the **Preview plots**.

## Add Plots to Workspace – Single Plot selection

- **Plots** are added sequentially to the currently displayed **Workspace** and their inserted position depends on the Workspace view mode **MLS** (“Workspace view modes” on page 126).
- **Plots** can be added to the **Workspace** multiple times.
- The indicator for plot addition persists with the **Workspace**.
- Single clicking a **Plot** adds it to the **Workspace** in the order described in “Create Workspace objects” on page 120.

## Add Plots to Workspace – Group selection

### Column header selection buttons

- Each column in the **Previews area** has an **Add column** button labeled with the parameter name positioned at the top of the column as a column header.
- By default, the parameter name is specified in the **Options dialog** (“Naming Options” on page 644).
- Clicking **Add column** adds the entire column of plots (excluding Histograms) to the **Workspace area** as described in “Create Workspace objects” on page 120.

### Row header selection buttons

- Each column in the **Previews area** has an **Add row** button labeled with the parameter name positioned at the beginning of the row as a row header.
- By default, the parameter name is specified in the **Options dialog** (“Naming Options” on page 644).
- Clicking **Add row** adds the entire row of plots (both Histograms and dual-parameter plots) to the **Workspace area** as described in “Create Workspace objects” on page 120.

### Select All split button

- Positioned at the top left corner of the **Previews area** is a split button, which can add the plots in either the upper right triangle or lower left triangle of the **Previews area** to the **Workspace**.
- Clicking the **top right** button adds the plots in the upper right triangle.
- Clicking the **lower left** button adds the plots in the lower right triangle.
- Plots are added to the **Workspace** as described in “Create Workspace objects” on page 120.

## Mouse over behavior

### Mouse over plot

- When the mouse pointer is hovered over a plot, a tooltip is displayed with the plot name and required action to add the plot to the Workspace.

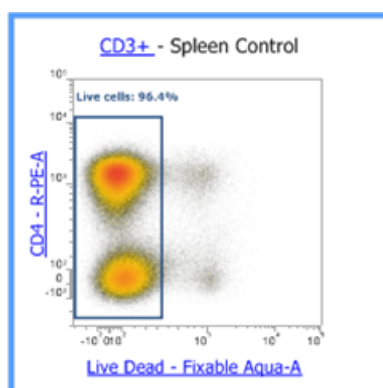
For single-parameter plots, the tooltip includes the X-parameter label.

FSC-A  
Click to add this plot to the workspace

For dual parameter plots the tooltip includes the X-parameter and Y-parameter labels in the format <Y Parameter> vs <X Parameter>.

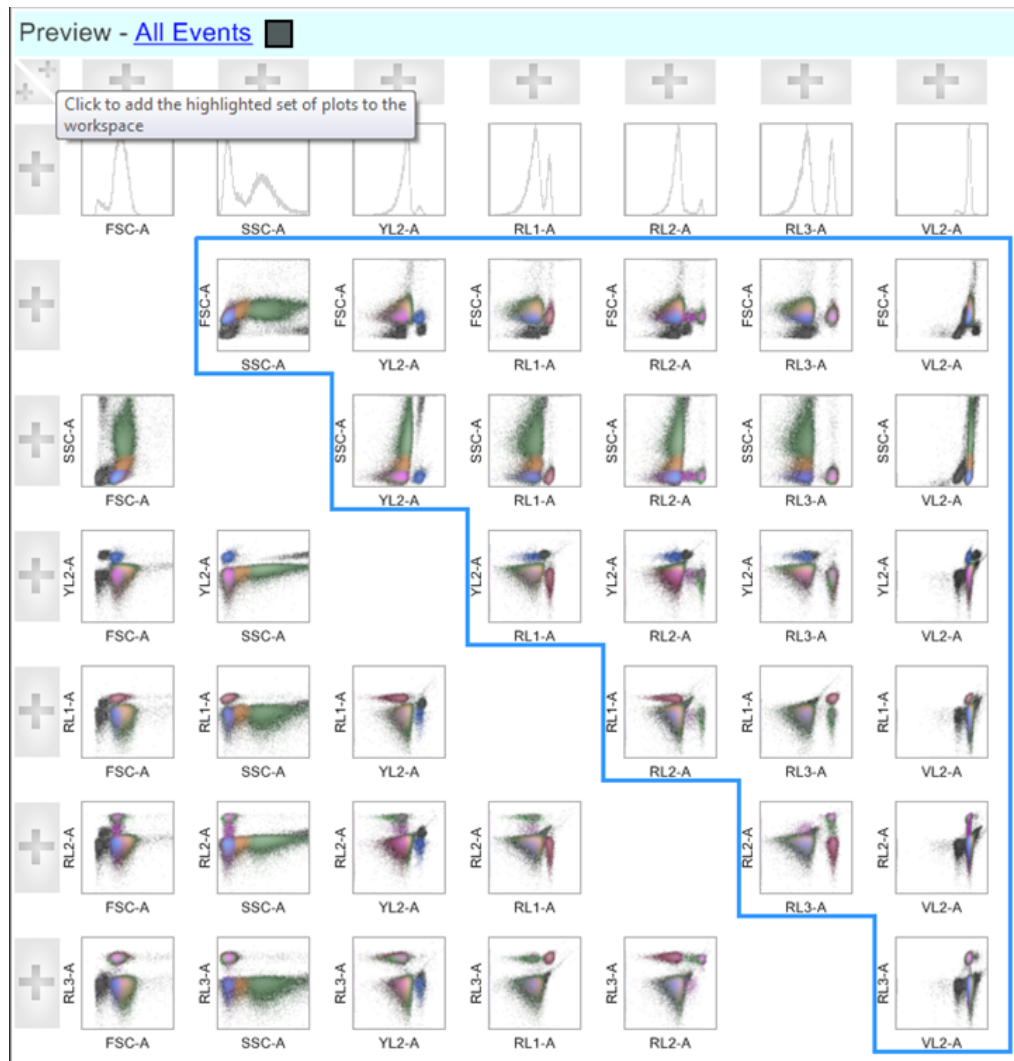
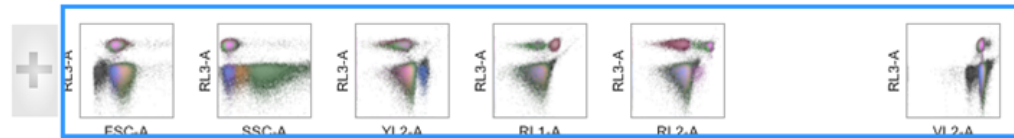
YL2-A vs FSC-A  
Click to add this plot to the workspace

- When the mouse pointer is hovered over a plot, a bounding rectangle is shown around the plot. (see “Select Workspace objects” on page 121).



## Mouse over buttons

- When the mouse pointer is hovered over **Group selection** buttons (i.e., column header selection, row header selection, or the **Select All** split buttons; “Add Plots to Workspace – Group selection” on page 135), the plots that will be added when the button is clicked are highlighted in an outline as described in “Workspace objects” on page 120.



- All group selection buttons show the following tooltip when the mouse pointer is hovered over them:

Click to add the highlighted set of plots to the workspace

# Plots

## Overview

The Attune™ Cytometric Software can show **Histogram**, **Dot**, **Density**, and **Precedence Density** plots on the **Workspace**.

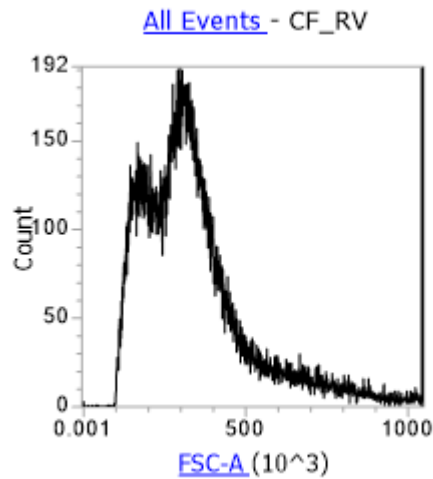
- Plots are created using the **Plots** group options present on the **Workspace** ribbon tab. The buttons in the **Plots** group are enabled when the **Workspace view** is in focus and at least two parameters are active, except for the **Histogram Plot** button, which requires only one active parameter.
- Plot options such as titles, axis labels, axis scaling, resolution, and colors used to show data points can be customized using the **Customize panel** (“Customize plot options” on page 421). Available customization options depend on the plot type selected.
- Default plot options are set using the **Options dialog** (“Fonts and Styles” on page 649).
- Dual-parameter plots (i.e., **Dot**, **Density**, and **Precedence Density** plots) show FSC values on the x-axis and SSC values on the y-axis by default.  
**Histogram** plots show FSC values on the x-axis by default.
- If FSC and/or SSC parameters are not selected for collection, the next parameter selected in the **Parameters** section of the **Instrument Settings panel** (“Parameters” on page 385) is used. The **Time** and **Event** parameters are only used if no other parameters are being collected.
- If the parameters that exist in the **Workspace** are disabled before recording data, a watermark on the plot states “Parameter(s) not available in dataset”. Any daughter plot also displays the watermark. The watermark does not appear until data are displayed or collected on the **Workspace**.
- Gates can be manipulated on top of the watermark. The plot scales are set to the last seen valid scale. Gate statistics from plots that do not have available parameters are displayed as **N/A**.



## Histogram plot

A **Histogram plot** is a graphical representation of single-parameter data and shows the relative number and distribution of events.

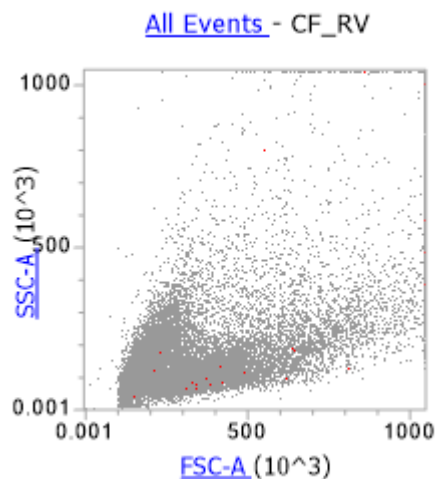
In a Histogram plot, the horizontal axis corresponds to the signal intensity of the selected parameter and the vertical axis represents the number of events (count).



- **Plot customization** options available for **Histograms plots** are **Resolution**, **Normalize Count**, **Use Shading**, and **Line Width**.
- The maximum plot resolution is 1024 channels, which is also the default resolution.
- The default line color for an ungated plot is set according to the selection in **Plot Options** in the **Options dialog** (“Plot Options” on page 653).
- The line color for a histogram plot that is based on a parent gate is determined by the parent gate’s color.

## Dot plot

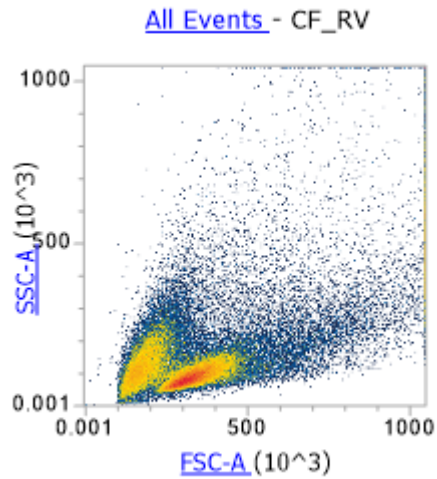
A **Dot plot** is a graphical representation of two-parameter data where each axis represents the signal intensity of one parameter. Each dot in the plot corresponds to one or more events detected above the threshold. Different colors are used to represent the parent gate of events that fall within bins on the plots.



- **Plot customization** options available for **Dot plots** are **Resolution** and **% of Events**.
- The maximum plot resolution is 1024 × 1024 channels. The default plot resolution is set at 256 × 256 channels.
- Where an event is in multiple gates, the precedence order determines which color is displayed. Where events from multiple gates fall within a single bin, the highest precedence gate color is displayed.
- Ungated events are painted grey by default. The default dot color for ungated events can be modified in **Plot Options** in the **Options dialog** (“Plot Options” on page 653).

## Density plot

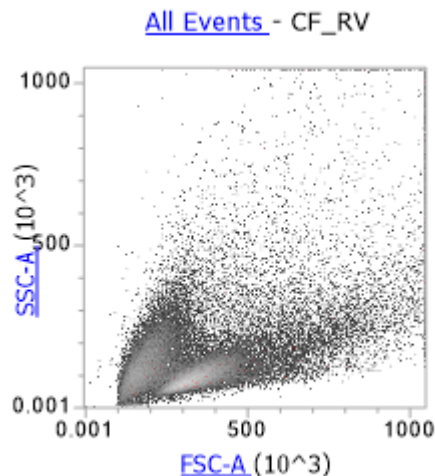
A **Density plot** is a graphical representation of two-parameter data where the colors represent the collection of events with the same intensity and each axis represents the signal intensity of one parameter.



- **Plot customization** options available for **Density plots** are **Resolution**, **Mode**, **Color**, and **% of Events**.
- The maximum plot resolution is 1024 × 1024 channels. The default plot resolution is set at 256 × 256 channels.

## Precedence density plot

A **Precedence Density plot** uses a combination of **Dot** and **Density** display, where a gradient indicates the number of events within each plot bin and color is used to show the parent gate of events present.



- **Plot customization** options available for **Precedence Density plots** are **Resolution**, **Mode**, and **% of Events**.
- The maximum dot plot resolution is 1024 × 1024 channels. The default plot resolution is set at 256 × 256 channels.

## Adjust plots

When you select individual plots, more controls become available, depending on the selected plot and system status.

- When **Plot Compensation** button on the **Compensation ribbon tab** (“View tab” on page 77) is turned **ON**, you can adjust compensation directly on the selected plot (“On-Plot compensation adjustment” on page 602).
- If the selected plot has an axis where the scale is set to Hyperlog™, a **Hyperlog™ transition value slider** (🔲) becomes available on each axis that is set to Hyperlog™.

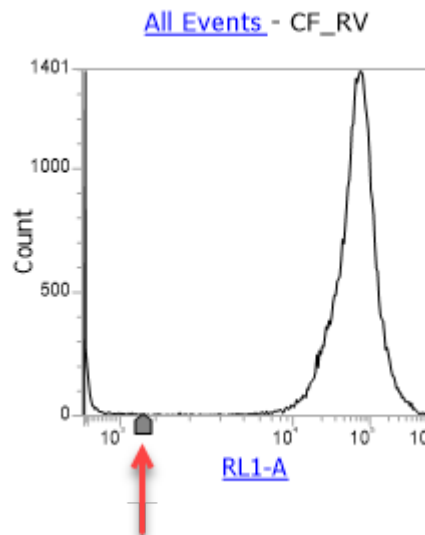


Figure 14 Histogram with Hyperlog™ slider

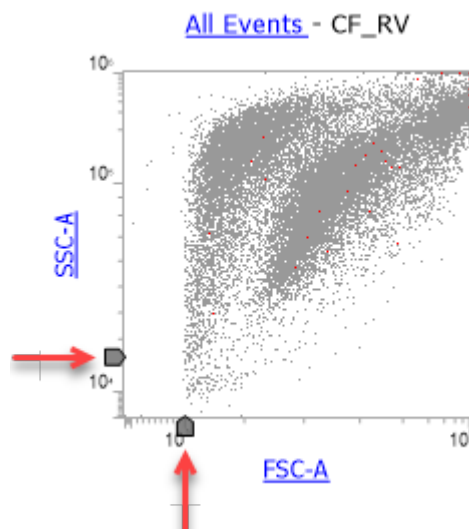


Figure 15 Dual-parameter plot with Hyperlog™ slider

- As you move the slider, the transform value is applied to the selected parameter on the selected plot (no other plots are updated). The value does not update in the customize panel when moving the slider.
- The slider position can be moved from 1 to 100,000.
- When you release the slider, the transition value is applied to all plots where the parameter is used, and the axis scaling changes to Hyperlog™.
- The new transform value is reflected in the **Plot Customize panel** for the selected plot and axis.
- The slider transform value applied to the data are linked to the axis value where the slider needle is depressed, and the axis tick marks are updated based on the newly applied transform.
- The slider needle is not displayed on plots when the workspace is printed, a plot is copied to the clipboard, or the plot is exported as an image.

# Gates

## Overview

**Gates** enable you to identify and analyze subsets of data (i.e., populations) by isolating a region in a selected plot. They can also be used to limit the number of events collected or stored.

After defining gates, you can combine them to create gates based on Boolean operators (i.e., Derived gates). You can use gated populations to generate statistics, show them in a hierarchal view, and create subsets within defined populations (i.e., daughter plots and associated statistics).

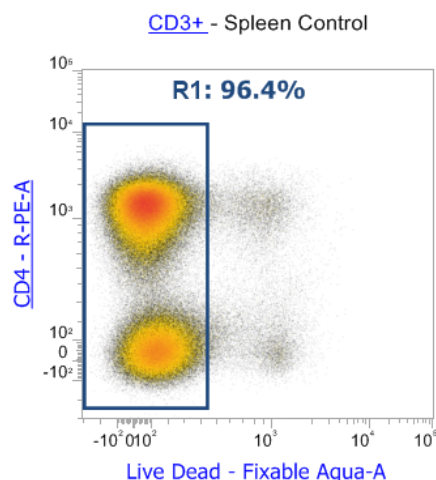
## Create and delete gates

- You can create gates on a selected plot using the **Gating Tools** available on the **Workspace ribbon tab**. Alternatively, you can use the **Plot context menu** (“Add gate” on page 192) to create gates.
- The following types of gates can be added on plots on a **Workspace**:
  - **Rectangular Gate** (“Rectangular Gate” on page 149)
  - **Oval Gate** (“Oval Gate” on page 150)
  - **Polygon Gate** (“Polygon Gate” on page 151)
  - **Quadrant Gate** (“Quadrant Gate” on page 152)
  - **Bent Quadrant Gate** (“Bent Quadrant Gate” on page 153)
  - **Histogram Gate** (“Histogram Gate” on page 155)
  - **Contour Autogate** (“Contour Autogate” on page 156)
  - **Ellipse Autogate** (“Ellipse autogate” on page 159)
  - **Derived Gate** (“Derived gate” on page 163)
- **Autogates** are defined by auto region bounding rectangles, which determine the autogate target area when the underlying data meet defined autogate criteria. If these criteria are met, an autogate of the specified type (contour or ellipse) is generated (see “Autogate gating rules” on page 162).
- Gate creation defaults, including autogates, are set in **Gate Options** in the **Options dialog** (“Gate Options” on page 656).
- You can create a maximum of 128 gates on plots on a **Workspace**, except for **Quadrant Gates**, of which at least 1000 can be created.
- You can create a maximum of 128 gating equations. A gating equation is generated when a new **Rectangle**, **Polygonal**, **Oval**, **Autogate**, or **Histogram Gate** is created, or when the **Derived Gate** option is used to create a compound gate equation.
- The maximum number of gates present in an individual gate equation is 28.
- When the **Workspace** reaches the maximum number of gates allowed, the relevant gate creation buttons on the **Workspace ribbon tab** and the options on the **Plot context menu** become disabled.
- To delete a gate, select it, and press the **Delete** key on the keyboard or the **Delete** button on the **Home tab**.

- When a gate is deleted, any **Derived Gate** that includes the deleted gate is modified to exclude the deleted gate.  
If the gate equation includes only the deleted gate, then the entire gate equation is deleted.
- Gate name, appearance, and position can be customized using the **Customize panel** (see “Customize gate options” on page 447).

## Gate name

- By default, each new gate is assigned a unique name, which consists of the letter R followed a number. The number increases incrementally for each new gate created, ensuring that each gate has a unique name (e.g., R1, R2,...).
- For **Quadrant gates**, the default gate name consists of the letter Q followed by an incrementally increasing number until a unique gate name is achieved (e.g., Q1, Q2,...).

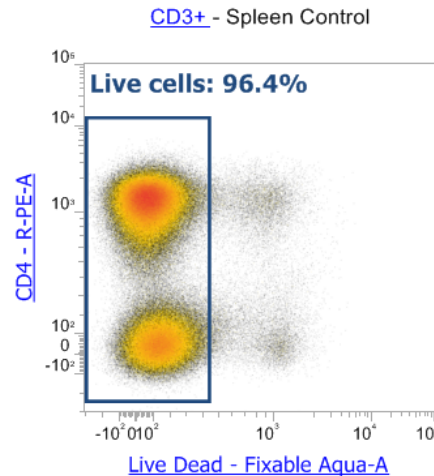


- The **gate name** and the associated **statistic** (optional) are displayed next to the gate on the relevant plot on separate lines. The font size auto-adjusts to optimally show the text.
- The display of a **statistic** is selected on the **Statistics ribbon tab** (“Statistics tab” on page 103) along with the desired statistic. Only one statistic can be shown and this applies to all plots. If a statistical value cannot be calculated, **N/A** is displayed instead of a numerical value.



- To change the assigned gate name, manually enter a name in the **Name** field in the **Customize panel** or select one of the predefined names from the **Quick Select** dropdown menu (“Name options” on page 459).

Alternatively, double-click the name of a selected gate to enter the name on plot edit mode.







- To reposition the gate name, first select the gate, then placing the cursor over the name. When the cursor changes to the standard precision select pointer (i.e., cross-hair), click and drag the gate name to a new position. You cannot drag the gate name outside the plot.
- You can move the gate name independently of the gate, but moving the gate keeps the gate name in the same relative position.

## Gate color

- When a gate is created, the **gate color** is automatically assigned.
- There are ten default colors defined for gates. As a new gate is created, the next available color is used. When an eleventh gate is created, the first color is reused. This process continues as each new gate is created.
- The **gate color** and the **opacity of the fill color** can be customized in the **Customize panel** (see “Customize gate options” on page 447).

## Move and resize gates

You can move and resize any gate. When you move or resize a gate, all populations derived from that gate are affected and statistics are automatically updated.

- When you click a gate to select it, the gate displays the **control points**, which are used for resizing or rotating the gate.
- When you move the mouse over the gate or the handle, the cursor changes its shape to indicate its action mode (**move**, **resize**, or **rotate**).
  -  **Move**: Used to move the selected object in any direction.
  -  **Resize**: Used to resize a single dimension (vertical or horizontal).
  -  **Diagonal resize**: Used to resize two dimensions simultaneously.
  -  **Rotate**: Used for freely rotating the selected object around a pivot point. **Rotate** function is available only for **Polygon** and **Rectangular gates**.
- To move a selected gate, grab it between the handles as indicated by the **Move** cursor, and then drag it to a new position. For opaque gates, you can also grab the center area of the gate to move it.
- You cannot move or resize **autogates** when they are **ON**. When the autogates are **OFF** or in the **ON – [NO FIT]** state, the underlying **bounding box** can be moved and resized.
- A **Quadrant gate** displays a **Move** cursor over the **intercept point**. For opaque quadrants, the center is not selectable for moving the gate.
- You cannot move a gate or quadrant intercept point outside of the plot axes, although a gate can exist outside the visible plot area, if an area that does not include the location of the gate was zoomed on the plot.

---

**Note:** It is also possible for a gate to exist outside of the plot limits, if the scale type or scale range of a plot has been modified after the gate has been defined (see “Off-plot gates” on page 164).

---

- When the **Rotate** icon is displayed on a selected gate, you can rotate the gate by dragging the indicated rotate control point.
- When a control point of a selected gate is being moved, the following **X** and **Y coordinates** are displayed in the **Application status bar**:
  - **Polygon Gate**: Coordinates of the point being moved are displayed.
  - **Histogram Gate**: Coordinates of the point being moved are displayed.
  - **Rectangular Gate**: Coordinates of the point being moved are displayed.  
If moving the top or right boundary line, the coordinates for the upper right control point are displayed.  
If moving the lower or left boundary line, the coordinates of the lower left control point are displayed.
  - **Oval Gate**: Coordinates of the center point are displayed.
  - **Autogates (Contour and Ellipse)**: Coordinates of the auto region bounding box are displayed.
  - **Quadrant Gate**: Coordinates of the intercept are displayed.
- When moving an entire **Rectangular**, **Polygon**, or **Oval gate**, or a **broken quadrant**, no coordinates are displayed.

## Rectangular Gate

To insert a **Rectangular gate** into a selected dual parameter plot, click the **Rectangular Gate** button on the **Workspace ribbon tab**.



Alternatively, select **Add Gate** ▶ **Rectangular** from the **Plot context menu**.

- To insert the gate, click the **Rectangular Gate** button to select, then click the desired dual parameter plot, which defines a corner of the rectangle.  
To draw the gate, drag out a rectangle from this point.
- When creating a Rectangular gate, the **Application status bar** displays the X and Y coordinates of the point being dragged.
- Select the **Rectangular Gate** button, then click a dual parameter plot to insert a default sized Rectangular gate centered on the point where the plot was clicked.  
The width of the Rectangular gate is 40% of the plot width and its default height is 20% of the plot width.
- On a single parameter plot, a no entry icon (⊘) indicates that a Rectangular gate cannot be created.
- **Rectangular gates** have eight **control points**.



- You can drag the center of each line forming the rectangle perpendicular to the line to resize the gate.
- You can drag the four corners of the rectangle, with the opposite corner being the anchored point. Its new position is determined by the rectangle formed by the anchored point and the final position where the mouse button is released.
- When you drag the rectangle, but not on a control point, the whole gate is dragged while maintaining its shape. For opaque gates, the center area is also selectable for moving the gate.

---

**Note:** If you are using an Attune™ CytPix™ Flow Cytometer, you can use **Rectangular gates** to set **Image gates**, which enable you filter the cell images displayed in **Image view** (see “Image capture gate” on page 276 for more information).

---

## Oval Gate

To insert an **Oval gate** into a selected dual parameter plot, click the **Oval Gate** button on the **Workspace ribbon tab**.

Alternatively, select **Add Gate** ▶ **Oval** from the **Plot context menu**.



- To insert the gate, click the **Oval Gate** button to select, then click the desired dual parameter plot, which defines a corner of the bounding box for the ellipse. To draw the gate, drag out an ellipse from this point.
- When an Oval gate is being created, the **Application status bar** displays the X and Y coordinates of the upper right corner of the bounding rectangle.
- Select the **Oval Gate** button, then click a dual parameter plot to insert a default sized Oval gate centered on the point where the plot was clicked. The default width of the Oval gate is 40% of the plot width and its default height is 20% of the plot height. The default Oval gate is created with the major axis horizontal.
- On a single parameter plot, a no entry icon (⊘) indicates that an Oval gate cannot be created.
- **Oval gates** have four **control points**. These represent the major and minor axis of the ellipse.
- Drag the control point in the direction of an axis to resize that axis.



- When you drag on the ellipse, but not on a control point, the whole gate is dragged while maintaining its shape. For opaque gates, the center area is also selectable for moving the gate.

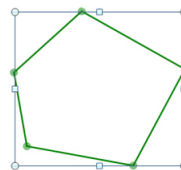
## Polygon Gate

To insert a **Polygon gate** into a selected dual parameter plot, click the **Polygon Gate** button on the **Workspace ribbon tab**.



Alternatively, select **Add Gate** ▶ **Polygon** from the **Plot context menu**.

- To insert the gate, click the **Polygon Gate** button to select, then click a dual parameter plot, which determines where the first point of the gate is placed. Continue to click the plot to add more points to the polygon to draw the gate around the population of interest. To complete the process, click the original point or double-click when inserting the last point.
- If there are at least three points, you can also complete the polygon by right clicking, which adds a line from the last to first point.  
If you right click when only two points have been defined, the gate is removed from the plot.
- The maximum allowable number of defined points on a Polygon gate is 40.
- When setting the first point of the polygon, a no-entry icon (⊘) is displayed when hovering outside a plot window. On a single parameter plot, a no entry icon indicates that a Polygon gate cannot be created.
- When a Polygon gate is being drawn, the **Application status bar** displays the X and Y coordinates of the cursor as each point is created.
- When creating a point of the polygon (other than the first point), if you move the cursor outside of the plot but remain within the application window (or child windows), a **precision select pointer** (i.e., **cross-hair**) continues to be displayed. If you click while the cursor is outside the plot, a control point is created at the coordinate closest to the clicked point.
- If you click outside the application window (or application child windows), the current gate being created is removed and the create mode is canceled.
- Clicking when the no-entry icon is displayed to create the first point of the polygon cancels the create mode without creating the point.
- Each vertex of the polygon is a control point. A bounding rectangle is displayed when the completed Polygon gate is selected.

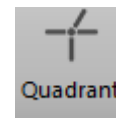


- Dragging a vertex moves that vertex only within the polygon.
- Dragging on a line within the polygon or the bounding rectangle moves the gate while maintaining the shape of the gate.
- Double-click a line within the polygon to insert a new vertex.
- Drag the line that forms the bounding rectangle perpendicular to the line to resize the gate.
- You can drag the four corners of the bounding rectangle, with the opposite corner being the anchored point.

- When you drag the bounding rectangle, but not on a vertex control point, or when you drag the line between vertex control points, the whole gate is dragged while maintaining its shape. For opaque gates, the center of the Polygon gate is also selectable for moving the gate.
- The default Polygon gate is defined by a bounding rectangle with 5 control points in the center of the plot. The width of the default bounding rectangle is 40% of the plot width and its default height is 20% of the plot height.

## Quadrant Gate

To insert a **Quadrant gate** into a selected dual parameter plot, click the **Quadrant Gate** button on the **Workspace ribbon tab**.



Alternatively, select **Add Gate** ▶ **Quadrant** from the **Plot context menu**.

- Only a single Quadrant gate can be created on a plot.
- Quadrant gates are formed from two **divider lines** (one vertical, one horizontal). To insert the gate, click the **Quadrant Gate** button, then click a point in the desired dual parameter plot. The point that is clicked in the plot determines where the **quadrant intercept point** is placed.
- On a single parameter plot, a no entry icon (⊘) indicates that a Quadrant gate cannot be created.
- After you have created the gate, you can edit the placement of the gate using the **control point** displayed on the **quadrant intercept**.
- By default, quadrants are added to a plot with the center point at the location that the plot is clicked.
- The quadrant is added orthogonally.
- By default, **quadrants** have five **control points**; one situated at the **divider intercept point** and one situated at each end of the two **dividers**.

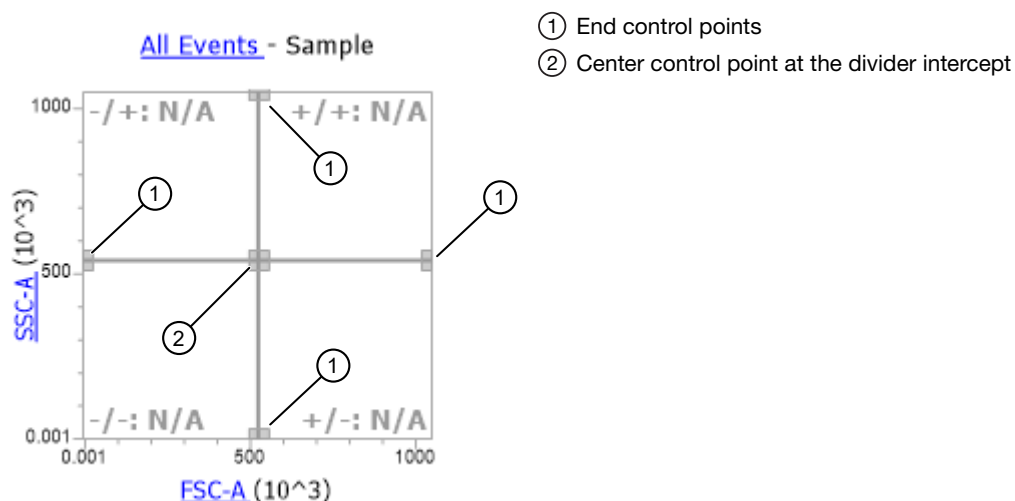


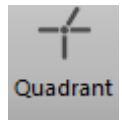
Figure 16 Quadrant gate control points

- While a **control point** is being moved, the X and Y coordinates of the intercept point are displayed in the **Application Status bar**.
- The quadrant intercept cannot be moved outside of the plot axes.

- The quadrant remains orthogonal until an **end control point** is moved (or the **Bent Quad** mode is selected in the **Quadrant gate customize panel**).
- If an end point is not moved, moving the center point moves the quadrant while maintaining the orthogonal nature of the quadrant gate.
- When an end point is moved, the quadrant becomes bendy and the **Bent Quad** button is selected in the **Customize Gate panel**.
- Any time a **quad vertex** is clicked with the left mouse button, the coordinates are displayed in the **Application Status bar**.

## Bent Quadrant Gate

To insert a **Bent Quadrant gate** into a selected dual parameter plot, click the **Quadrant Gate** button on the **Workspace ribbon tab**, then insert a **Quadrant gate** in the plot. After the **Quadrant gate** is inserted, move an **end control point** to make the gate a **Bent Quadrant gate** (or select the **Bent Quadrant** option in the **Customize Gate panel**).



Alternatively, select **Add Gate** ▶ **Quadrant** from the **Plot context menu** (“Insert plot statistics” on page 192), then make the gate a **Bent Quadrant gate** as described above.

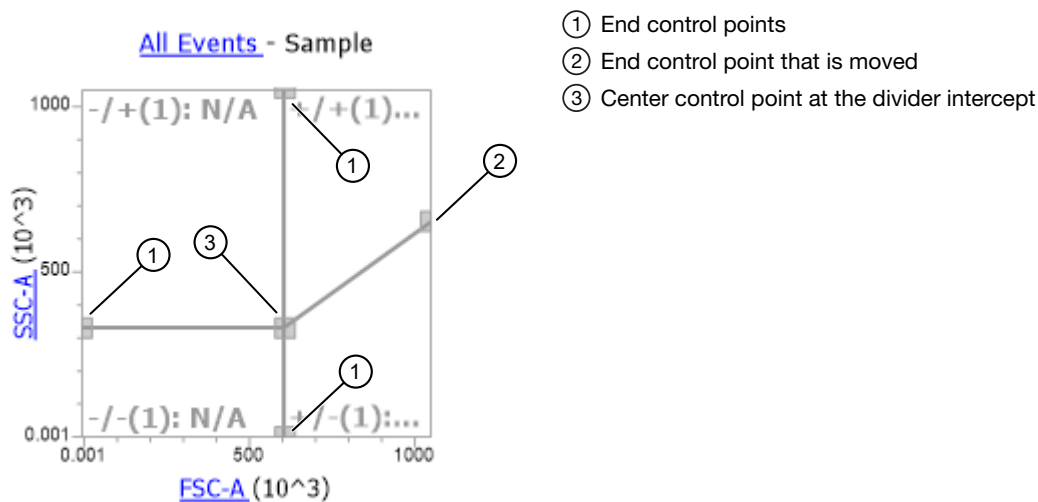


Figure 17 Quadrant gate control points. Moving an end control point turns the Quadrant gate into a Bent Quadrant gate.

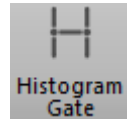
- In a **Bent Quadrant gate**, moving the **center point** only moves the center point while the **end points** remain at their original positions.
- The coordinates of the end points of a quadrant are extrapolated from the center point to a limit of PnR for positive coordinates and  $-PnR$  for negative coordinates. You can only move an end point within these limits.
- You cannot move an end point from its corresponding border (i.e., the left end point remains on the left boundary of the plot, the top on the top, the right on the right, and the bottom on the bottom) where the coordinate of that border is fixed at its limits (for example, the left end point has an x-coordinate of  $-PnR$ ).

- For a **Bent Quadrant gate**, the status bar displays the following coordinates:
  - **Top X**: User defined within plot limits for X axis
  - **Top Y**: Plot's Y Max
  - **Right**: Plot's X Max
  - **Right Y**: User defined within plot limits for Y axis
  - **Left X**: Plot's X Min
  - **Left Y**: User defined within plot limits for Y axis
  - **Bottom X**: User defined within plot limits for X axis
  - **Bottom Y**: Plot's Y Min
  - **Center Points**: User defined within plot scale space (visible and not visible) (i.e., **Linear**, **Log**, or **Hyperlog™**)
- When the plot scale and range is changed where the center point is outside the visible bounds of the plot, the quadrant is displayed in red to indicate that the quadrant center point is outside the visible plot space. If the range and scale are changed so that the center point is visible, it is rendered without red and with the center point in its specified location.
- If the range or scale is changed such that a quadrant vertex is no longer on its corresponding plot border, it is drawn with a dashed line.



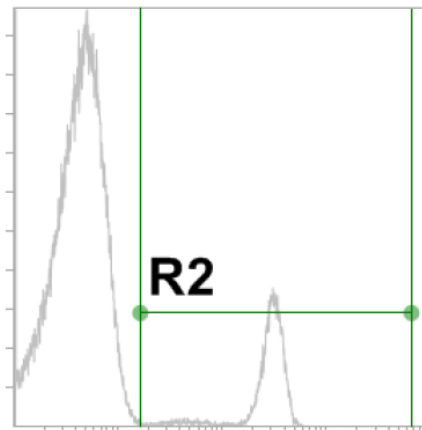
## Histogram Gate

To insert a **Histogram gate** into a selected single parameter plot, click the **Histogram Gate** button on the **Workspace ribbon tab**.



Alternatively, select **Add Gate** ▶ **Histogram** from the **Plot context menu**.

- To insert the gate, click the **Histogram Gate** button, then click the desired single parameter plot and drag out a line to determine the full gate size.
- When a Histogram gate is being created, the **Application status bar** displays the X and Y coordinates of the point being dragged.
- Selecting the **Histogram Gate** button and clicking on a single parameter plot inserts a default sized Histogram gate centered on the point where the plot was clicked. The default width of the Histogram gate is 40% of the plot width. Full height vertical lines are used to mark the upper and lower limits of the gate.
- On a dual parameter plot, a no entry icon (⊘) indicates that a Histogram gate cannot be created.
- Histogram gates have **control points** situated at each end of the gate. When the gate is selected, **control points** are displayed on each boundary.



- Dragging the control points or the vertical boundary line moves the end points in the direction of the appropriate axis. Dragging the horizontal line, but not on a control point, enables the whole gate to be moved.
- The Histogram gate cannot be moved outside the plot axes. For opaque gates, the center is also selectable for moving the gate.
- When needed, the default position of the Histogram gate is the center point of the gate set to the center of the plot.

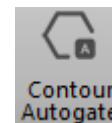
---

**Note:** If you are using an Attune™ CytPix™ Flow Cytometer, you can use **Histogram gates** to set **Image gates**, which enable you filter the cell images displayed in **Image view** (see “Image capture gate” on page 276 for more information).

---

## Contour Autogate

To insert a **Contour autogate** into a selected dual-parameter plot, click the **Contour Autogate** button on the **Workspace ribbon tab**.



This enables you to draw a rectangular region (**auto region bounding rectangle**), which becomes the basis of the **Contour autogate target area**.

Alternatively, select **Add Gate** ▶ **Contour Autogate** from the **Plot context menu**.

---

**Note:** **Autogates** are set if the criteria for population identification are met within a defined **auto region bounding rectangle** (see “Autogate gating rules” on page 162).

---

- To insert the **Contour auto region bounding rectangle**, click the **Contour Autogate** button, then click the desired dual parameter plot and drag out a line to determine the full gate size.
- When a Contour autogate is being created, the **Application status bar** displays the X and Y coordinates of the point being dragged.
- Selecting the **Contour Autogate** button and clicking on a plot inserts a default sized **Contour auto region bounding rectangle** that is centered on the point where the plot was clicked. The default width and height of an autogating region is 20% of the plot width and height.
- A no entry icon (⊘) indicates that a Contour autogate cannot be created in that area of the Plots view.
- When the Contour auto bounding region is first created, it is in the **unautogated** state, which is indicated by dashed lines that make up the rectangle (see “Contour auto region states” on page 157).

If the autogating criteria are met when the gate is **ON**, then the auto region bounding rectangle automatically forms an autogate.

- When a successful autogate forms in the **ON** state, a smaller bounding rectangle that reflects the new Contour autogate is created, and the Application status bar displays the X and Y coordinates of the new bounding rectangle.

However, the coordinates of the bounding box for an autogate that is in the **ON – [NO FIT]** or the **OFF** state reflect the original, larger auto region bounding box.

- Contour auto region bounding rectangles have eight control points.
- If the autogate is **OFF** or in the **ON – [NO FIT]** state, you can drag the center of each line forming the rectangle perpendicular to the line to resize the auto region.

If the autogate is in the **ON** state, you cannot resize the auto region.



- You can drag the four corners of the rectangle, with the opposite corner being the anchored point. Its new position is determined by the rectangle formed by the anchored point and the final position where the mouse button is released.

- When you drag the rectangle, but not on a control point, the whole auto region is dragged while maintaining its shape. For opaque gates, the center area is also selectable for moving the gate.
- To delete a gate, select it, then press the **Delete** key or select **Delete** in the right-click menu after the gate is selected.

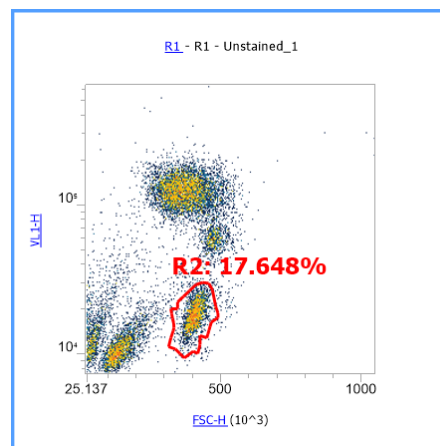
**Note:** Auto region default settings are dictated by the setting in the **Gate options** (“Gate Options” on page 656).

## Contour auto region states

**State:** Autogate **ON**

**Visualisation:** Solid line smoothed polygon

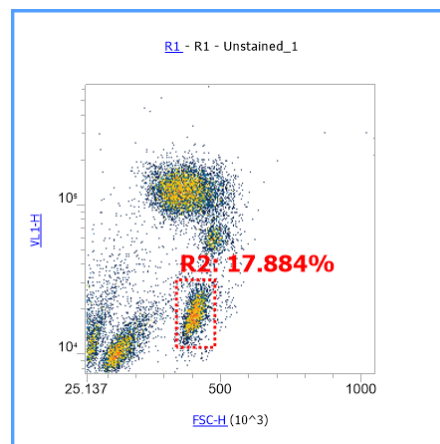
**Manipulation options:** Control points displayed when selected. Size and position are not manipulatable.



**State:** Autogate **OFF**

**Visualisation:** Dashed line polygon

**Manipulation options:** Control points displayed when selected. Size and position are manipulatable like a rectangular gate.



**State:** Autogate **ON** – [NO FIT]

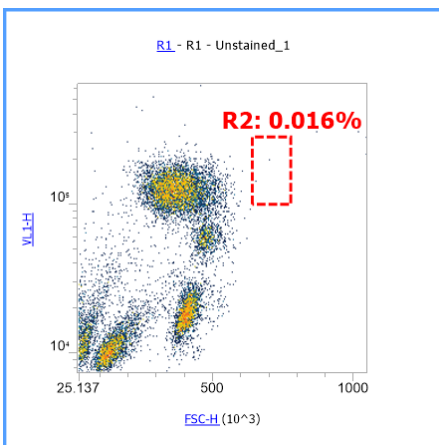
**Visualisation:** Dashed line rectangle

**Manipulation options:** Control points displayed when selected. Size and position are manipulatable based on rectangular region rules.

---

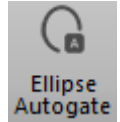
**Note:** After manipulation, the Auto Region attempts to autogate in this state. If successful, its state changes to **ON**.

---



## Ellipse autogate

To insert an **Ellipse autogate** into a selected dual-parameter plot, click the **Ellipse Autogate** button on the **Workspace ribbon tab**.



This enables you to draw a rectangular region (**auto region bounding rectangle**), which becomes the basis of the **Ellipse autogate target area**.

Alternatively, select **Add Gate** ▶ **Ellipse Autogate** from the **Plot context menu**.

---

**Note:** **Autogates** are set if the criteria for population identification are met within a defined **auto region bounding rectangle** (see “Autogate gating rules” on page 162).

---

- To insert the **Ellipse auto region bounding rectangle**, click the **Ellipse Autogate** button, then click the desired dual parameter plot and drag out a line to determine the full gate size.
- When an Ellipse Autogate is being created, the **Application status bar** displays the X and Y coordinates of the point being dragged.
- Selecting the **Ellipse Autogate** button and clicking on a plot inserts a default sized **Ellipse auto region bounding rectangle** that is centered on the point where the plot was clicked. The default width and height of an autogating region is 20% of the plot width and height.
- A no entry icon (⊘) indicates that an Ellipse autogate cannot be created in that area of the Plots view.
- When the Ellipse auto bounding region is first created, it is in the **unautogated** state, which is indicated by dashed lines that make up the rectangle (see “Ellipse auto region states” on page 160).

If the autogating criteria are met when the gate is **ON**, then the auto region bounding rectangle automatically forms an autogate.

- When a successful autogate forms in the **ON** state, a smaller bounding rectangle that reflects the new Ellipse autogate is created, and the Application status bar displays the X and Y coordinates of the new bounding rectangle.

However, the coordinates of the bounding box for an autogate that is in the **ON – [NO FIT]** or the **OFF** state reflect the original, larger auto region bounding box.

- Ellipse auto region bounding rectangles have eight control points.
- If the autogate is **OFF** or in the **ON – [NO FIT]** state, you can drag the center of each line forming the rectangle perpendicular to the line to resize the auto region.



If the autogate is in the **ON** state, you cannot resize the auto region.

- You can drag the four corners of the rectangle, with the opposite corner being the anchored point. Its new position is determined by the rectangle formed by the anchored point and the final position where the mouse button is released.

- When you drag the rectangle, but not on a control point, the whole auto region is dragged while maintaining its shape. For opaque gates, the center area is also selectable for moving the gate.
- To delete a gate, select it, then press the **Delete** key or select **Delete** in the right-click menu after the gate is selected.

---

**Note:** Auto region default settings are dictated by the setting in the **Gate options** (“Gate Options” on page 656).

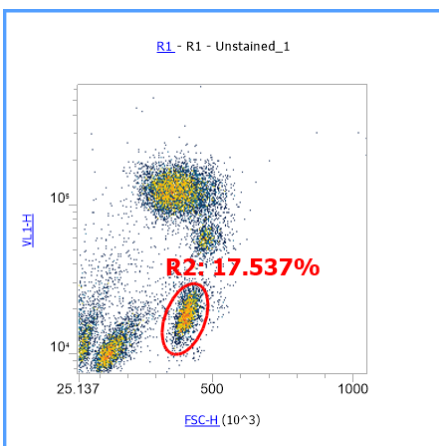
---

## Ellipse auto region states

**State:** Autogate **ON**

**Visualisation:** Solid line smoothed ellipse

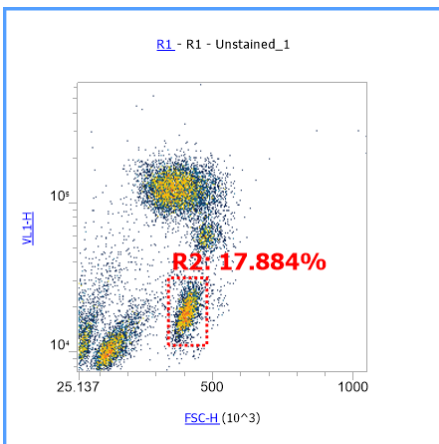
**Manipulation options:** Control points displayed when selected. Size and position are not manipulatable.



**State:** Autogate **OFF**

**Visualisation:** Dashed line ellipse

**Manipulation options:** Control points displayed when selected. Size and position are manipulatable like a rectangular gate.



**State:** Autogate **ON** – **[NO FIT]**

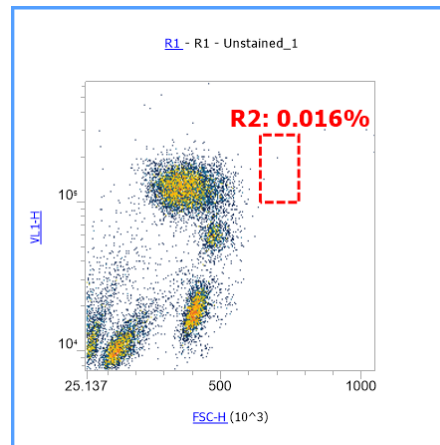
**Visualisation:** Dashed line rectangle

**Manipulation options:** Control points displayed when selected. Size and position are manipulatable based on rectangular region rules.

---

**Note:** After manipulation, the Auto Region attempts to autogate in this state. If successful, its state changes to **ON**.

---



## Autogate gating rules

Autogates are defined by **auto region bounding rectangles**, which determine the **autogate target area** for underlying data based on autogate settings. If the criteria defined in autogate settings are met, the auto region of the specified type (**contour** or **ellipse**) is created in the bounding rectangle.

The following table describes the potential auto region states at any given time after creation.

State	Description
Autogate <b>ON</b>	Auto region has been successfully autogated around a population or populations. The auto region is represented as the appropriate autogate type ( <b>Ellipse</b> or <b>Contour</b> ).  <b>ON</b> indicates that the auto region will attempt to autogate under any of the reautogate scenarios described in this section.
Autogate <b>ON – [NO FIT]</b>	Auto region has not been successfully autogated. The auto region is represented as a rectangular region type with a dashed boundary line.  <b>ON</b> indicates that the auto region will attempt to autogate under any of the reautogate scenarios described in this section.
Autogate <b>OFF</b>	Auto region has not been successfully autogated. The auto region is represented as a rectangular region with a dotted boundary line.  <b>OFF</b> indicates the auto region will <b>not</b> attempt to autogate under any of the reautogate scenarios described in this section.

### Re-autogating

An auto region with the Autogate ON setting will re-attempt to autogate based on its auto region bounding rectangle co-ordinates and its current autogate settings when new data are acquired, when FCS files are imported in a shared workspace, or when present in an imported workspace and applied to underlying data.



## Derived gate

**Derived gates** are gates based on **Boolean operators** applied to existing gates. You can create a derived gate when there is at least one gate present in the **Workspace**, with any or no plot selected.

- You can create a **Derived gate** using the **Derived Gate dialog** (“Derived gate dialog” on page 739).

To open the **Derived Gate dialog**, click the **Derived Gate** button on the **Workspace ribbon tab**.



Alternatively, select **Add Gate** ▶ **Derived Gate** from the Plot context menu.

- The **Boolean operators** available for creating Derived gates are **blank**, **AND**, **OR**, **XOR**, **AND NOT**, and **OR NOT**.

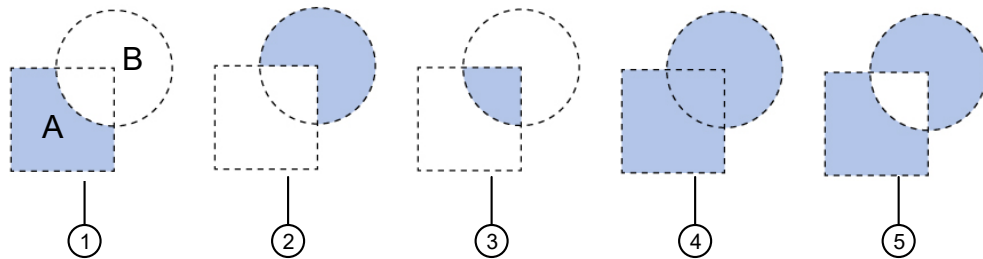


Figure 18 Boolean operators available for creating Derived gates.

- ① A NOT B
  - ② B NOT A
  - ③ A AND B (intersected)
  - ④ A OR B (joined)
  - ⑤ A XOR B
- For more information about creating **Derived gates**, see “Derived gate dialog” on page 739.

## Off-plot gates

It is possible for a gate to exist outside of the plot limits, if the scale type (linear, logarithmic, or Hyperlog™) or range (minimum and maximum of the plot axes) of a plot is modified after the gate has been defined. Gates are termed "off-plot", if their coordinates do not exist within the scale type or range. Off-plot gates can be partially or fully off-plot.

- If the gate is partially off the visible plot scale range and the gate coordinates exist within the plot scale type, the displayed gate is clipped at the plot axis or plot boundary.

① Gates partially off the plot

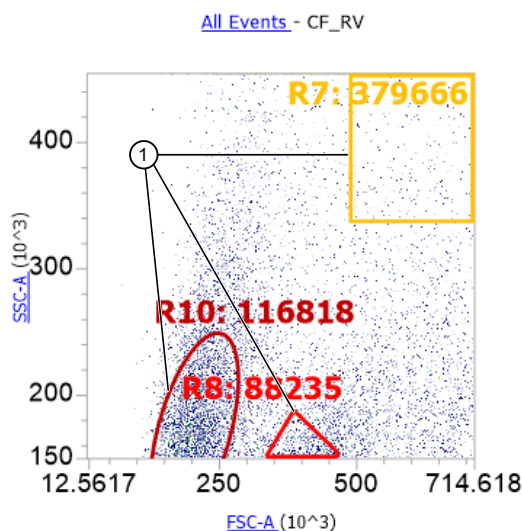


Figure 19 Gates partially off the plot boundary

- If the gate is partially off the visible plot scale range and the gate coordinates do not exist within the plot scale type (for example, after changing from Linear scale to Log scale with negative coordinates), the gate is shown with the control points depicting the undefined coordinates filled with red, when selected.

① Gate coordinates are undefined in Log scale – the control points are filled with red.

② Gate coordinates are within the plot scale range – the control points are filled with white.

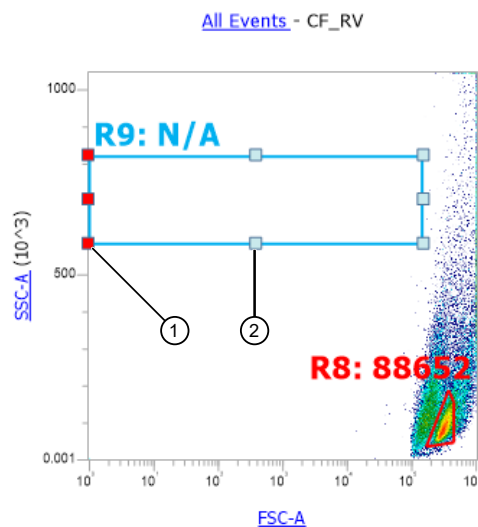
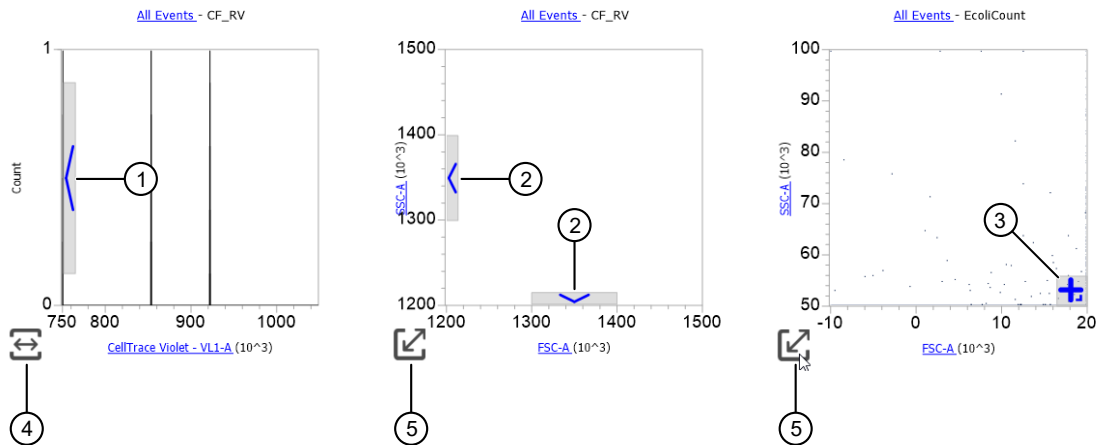


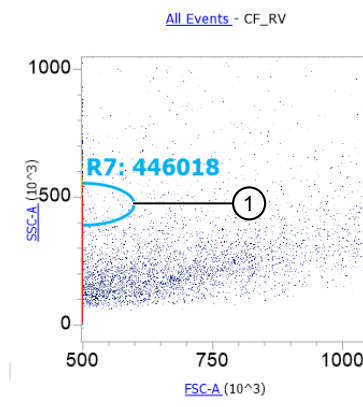
Figure 20 Gate with undefined gate coordinates

- Dragging the gate away from plot axis moves the gate into the defined space.
- For a **Histogram gate** on a Histogram plot, and **Rectangular, Polygon, Autogates, and Quadrant gates** on two-parameter plots, if one or more gates are completely off-plot, the plot shows one or more **off-plot gate indicators** (depending on the number and position of the off-plot gates) and an **axis expansion button**, so that the presence and location of the off-plot gate can be ascertained.



**Figure 21** Gate completely off-plot – Histogram gate (left) on a Histogram plot, Rectangular, Polygon, or Autogate on a two-parameter plot (middle), and Quadrant gate on a two-parameter plot (right).

- ① Off-plot gate indicator for a Histogram gate. The gate is off-plot to the left.
  - ② Off-plot gate indicators for Rectangular, Polygon, or Autogates on two-parameter plots. The gates are off-plot to the left and bottom of the plot.
  - ③ Off-plot gate indicator for a Quadrant gate on a two-parameter plot. The quadrant center falls off-plot to the lower right of the plot.
  - ④ Axis expansion button for Histogram plots.
  - ⑤ Axis expansion button for two-parameter plots.
- Plots with off-plot **Oval gates** do not show off-plot gate indicators. Instead, the gate is partially displayed next to the plot axis for the off-plot parameter.



**Figure 22** Gate completely off the plot – Oval gate

- ① Oval gate completely off plot – gate is displayed next to the plot axis with the default size for the off-plot parameter.

### Off-plot gate indicators - Histogram plots

If one or more histogram gates are completely off-plot, an **off-plot gate indicator** is displayed to the far left and/or the far right of the of the x-axis depending on the off-plot position of gate. In addition, an **axis expansion button** is displayed outside the plot axes at the left bottom of the plot.

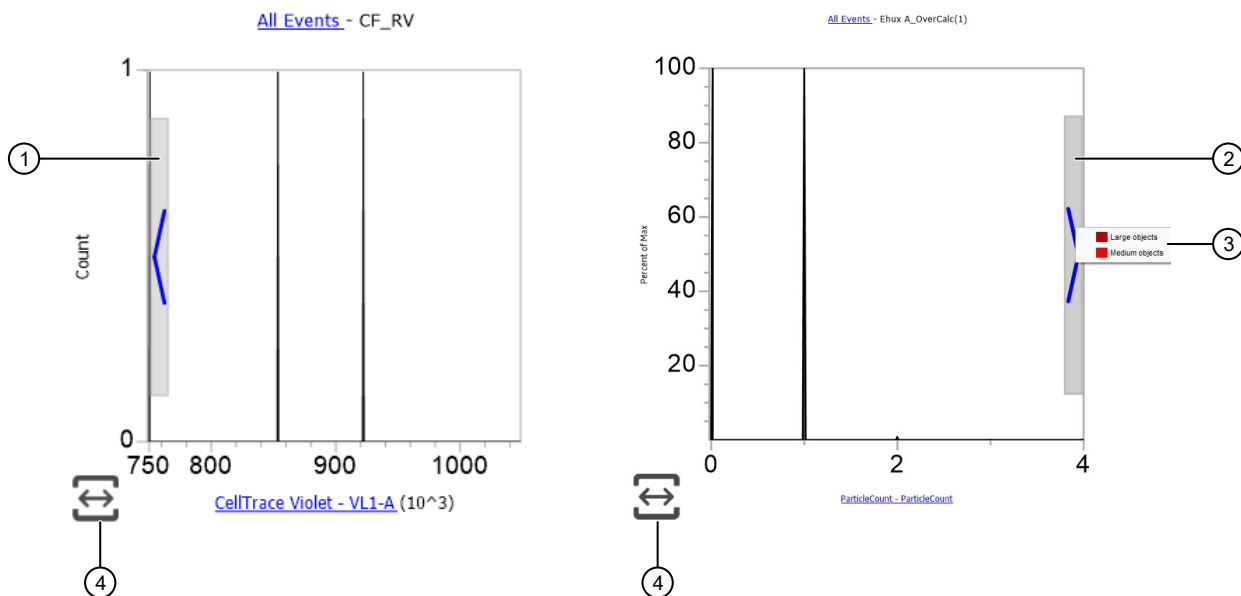


Figure 23 Off-plot gate indicators for Histogram plots

- ① Off-plot gate indicator (off-plot gate is to the left of the plot scale)
  - ② Off-plot gate indicator (off-plot gate is to the right of the plot scale)
  - ③ Off-plot gate indicator context menu
  - ④ Axis expansion button for Histogram plots
- Double-click the **axis expansion button** to expand the plot scale to the PnR/gate limit maximum.

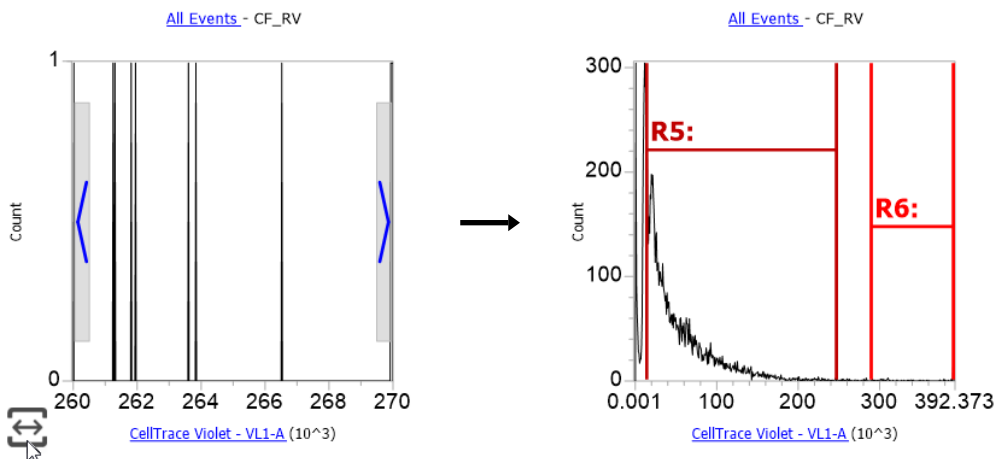


Figure 24 Use the axis expansion button to expand plot scale to include all off-plot gates.

- Click the **off-plot gate indicator** to expand the axis to the furthest boundary of the off-plot gate.

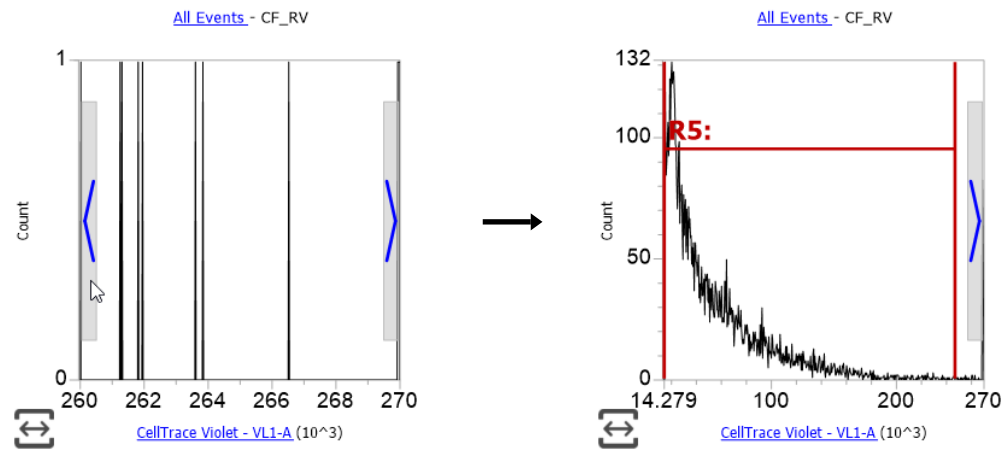


Figure 25 Use the off-plot gate indicator to expand the plot scale to include all off-plot gates in the indicator direction (in this example, the indicator that points to the left).

- Right-click the **off-plot gate indicator** to view the **off-plot gate indicator context menu**, which lists the off-plot gates alphabetically in the sort order R1, R2, R3 and so on. Select a gate from the context menu to expand the plot to the furthest boundary of the selected gate.

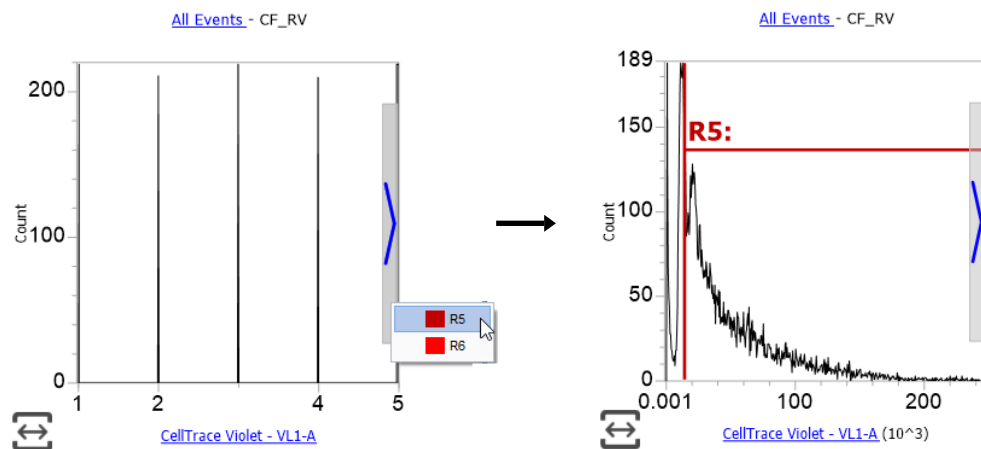


Figure 26 Use the off-plot gate indicator context menu to expand plot scale only to the boundary of the selected gate.

- When the **axis expansion button** or the **off-plot gate indicator** is clicked, the **Scale** mode is set to **Manual** and the plot scale is expanded to minimum and maximum values as determined by the gate/PnR bounds. This includes negative space for **Hyperlog™** and **Linear** scales.
- If a gate exists in the negative scale when a **Log** scale is used, the scale is changed to **Hyperlog™**. Any gates that are still off-plot continue to be represented by an **off-plot gate indicator**.
- Following plot expansion, if the axes are manually readjusted to cause one or more gates to show off-plot behavior, the **axis expansion button** and the **off-plot gate indicator** reappear on the plot.

## Off-plot gate indicators - Two-parameter plots (Rectangular, Polygon, and Autogates)

If one or more **Rectangular**, **Polygon**, or **Autogates** are completely off-plot, one or more **off-plot gate indicators** are displayed on the plot depending on the off-plot position of gates, with the **chevron** of the indicator pointing to the position of the off-plot gate. In addition, an **axis expansion button** is displayed outside the plot axes at the left bottom of the plot.

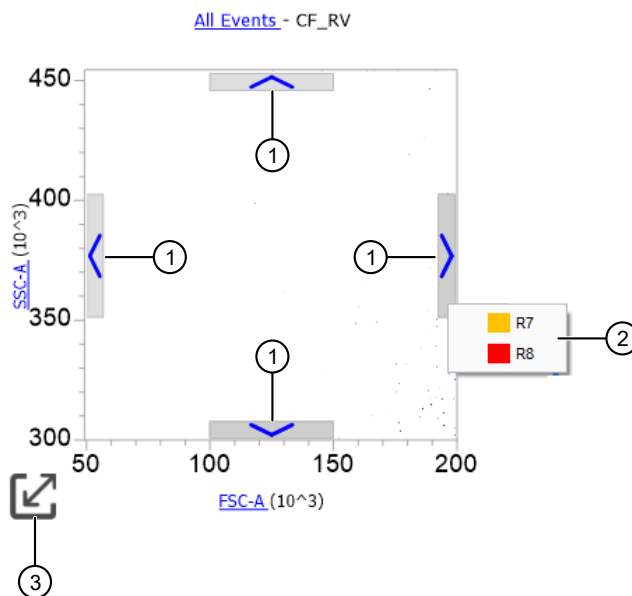


Figure 27 Off-plot gate indicators for two-parameter plots (Rectangular, Polygon, and Autogates)

- ① Off-plot gate indicator
- ② Off-plot gate indicator context menu
- ③ Axis expansion button for two-parameter plots
- Double-click the **axis expansion button** to expand the plot scale to the PnR/gate limit maximum.

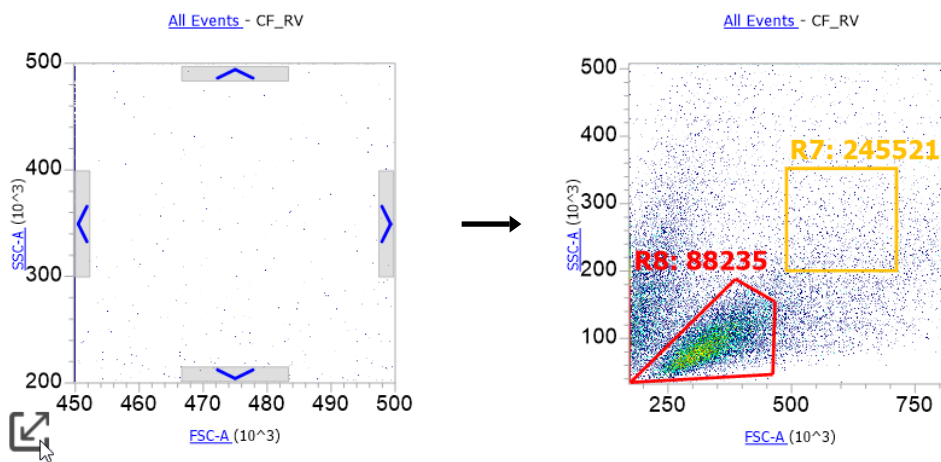


Figure 28 Use the axis expansion button to expand plot scale to include all off-plot gates.

- Click the **off-plot gate indicator** to expand the axis to the furthest boundary of the off-plot gate that is located in the direction as specified by the chevron of the indicator.

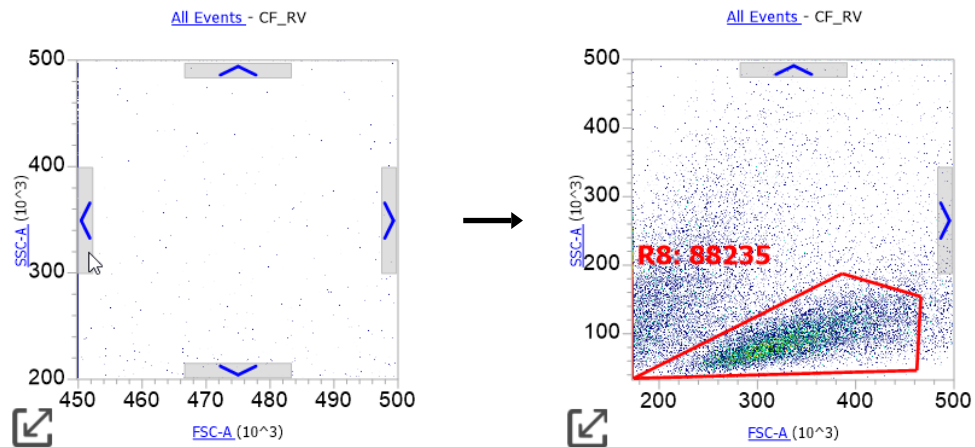


Figure 29 Use the off-plot gate indicator to expand the plot scale to include all off-plot gates in the indicator direction (in this example, the indicator that points to the left).

- Right-click the **off-plot gate indicator** to view the **off-plot gate indicator context menu**, which lists the off-plot gates alphabetically in the sort order R1, R2, R3 and so on. Select a gate from the context menu to expand the plot to the furthest boundary of the selected gate.

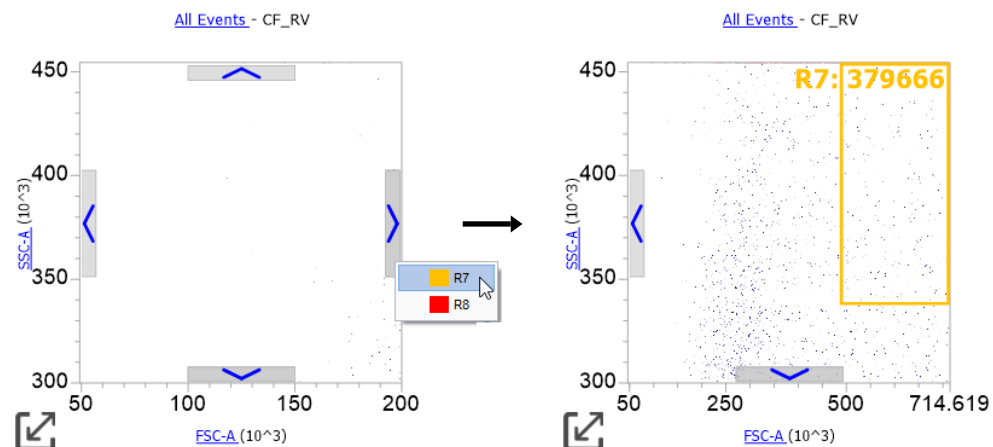


Figure 30 Use the off-plot gate indicator context menu to expand plot scale only to the boundary of the selected gate.

- When the **axis expansion button** or the **off-plot gate indicator** is clicked, the **Scale** mode is set to **Manual** and the plot scale is expanded to minimum and maximum values as determined by the gate/PnR bounds. This includes negative space for **Hyperlog™** and **Linear** scales.
- If a gate exists in the negative scale when a **Log** scale is used, the scale is changed to **Hyperlog™**. Any gates that are still off-plot continue to be represented by an **off-plot gate indicator**.
- Following plot expansion, if the axes are manually readjusted to cause one or more gates to show off-plot behavior, the **axis expansion button** and the **off-plot gate indicator** reappear on the plot.

## Off-plot gate indicators - Two-parameter plots (Quadrant gates)

If the center point of a **Quadrant gate** is off-plot, a cross-shaped **off-plot gate indicator** is displayed on the plot. In addition, an **axis expansion button** is displayed outside the plot axes at the left bottom of the plot.

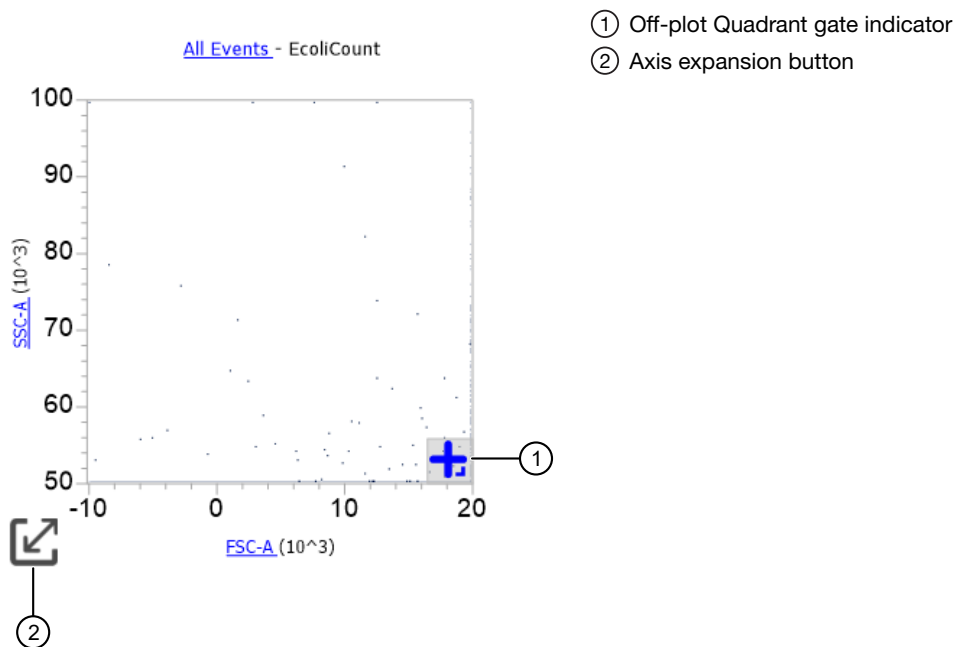


Figure 31 Off-plot gate indicators for Quadrant gates

- The position of the cross-shaped **off-plot gate indicator** reflects the location of the center of the off-plot quadrant gate and the **chevron** icon points to the relevant corner of the plot. For example, if the off-plot quadrant gate center falls into the top left section of the plot, the **off-plot gate indicator** is positioned in the top left of the plot. If the off-plot quadrant gate center falls into the top right section of the plot, the **off-plot gate indicator** is positioned in the top right of the plot, and so on. If the off-plot quadrant gate center falls directly on one of the intersect lines, the **off-plot gate indicator** is shown in both sections. For example, if the off-plot quadrant center is off plot and aligned with the left intersect, the **off-plot gate indicator** is shown in the top and lower left corners of the plot. Both indicators will have identical behavior for axes expansion.



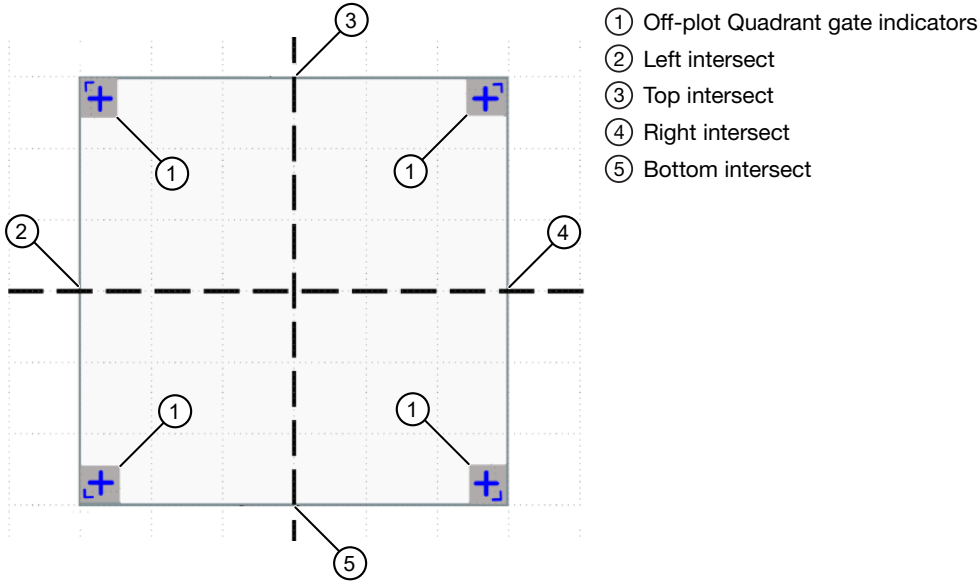


Figure 32 Off-plot Quadrant gate indicator behavior

- Double-click the **axis expansion button** to expand the plot scale to the PnR/gate limit maximum.

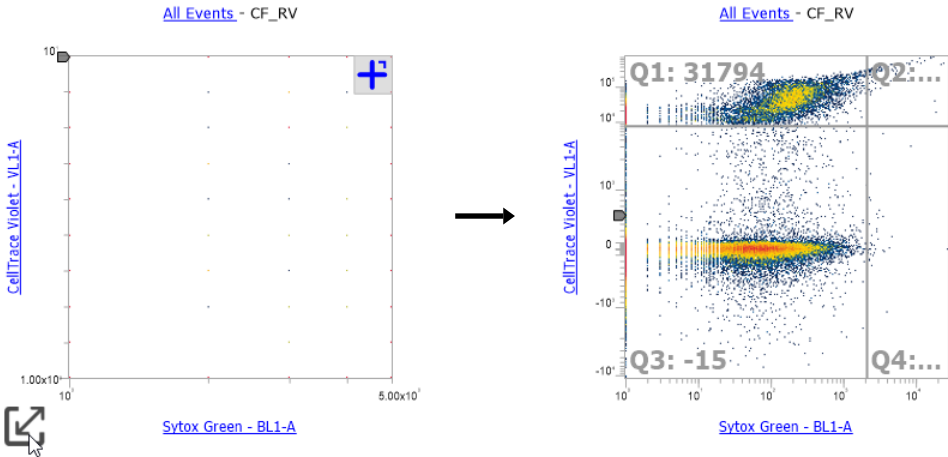


Figure 33 Use the axis expansion button to expand plot scale to include the off-plot Quadrant gate.

- Click the **off-plot gate indicator** to expand the plot axes to encompass PnR maximum and position the center of the quadrant gate in the middle of the plot (if possible).

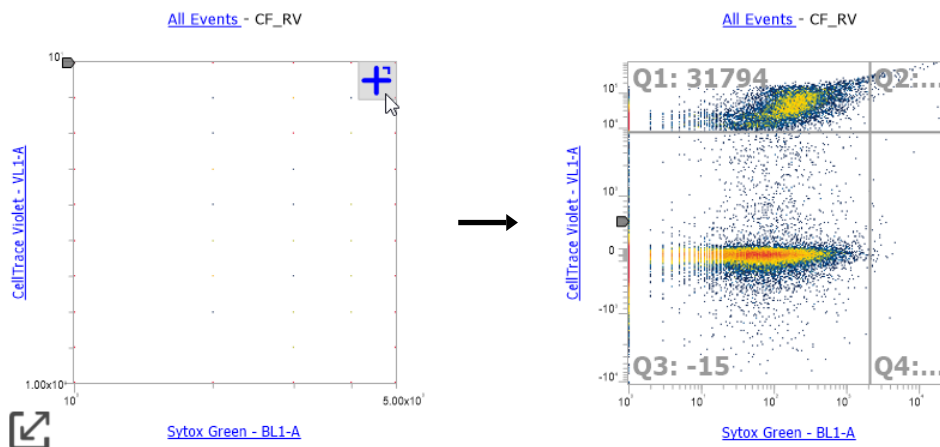


Figure 34 Use the off-plot gate indicator to expand the plot scale to include the off-plot Quadrant gate.

- Following plot expansion, if the axes are manually readjusted to cause the gate to show off-plot behavior, the **axis expansion button** and the **off-plot gate indicator** reappear on the plot.

## Gate behavior

When the **plot parameter number** or **plot parameter range** is changed, the gate is resized or deleted according to the rules described below.

When the **plot scale range**, **scale type**, **plot axis parameters**, or **plot type** is changed, the gates are resized or deleted as described below.

### Changing plot scale range

- If the **plot scale range** is changed on any axis, all gates remain fixed in their specified locations.
- If the change in scale range results in a gate being off plot, the gate behaves as described in “Off-plot gates” on page 164.

### Changing plot scale type

- If the **plot scale type** is changed (**Linear**, **Log**, **Hyperlog™**) on any axis, the coordinates of the vertices for all gates (excluding Oval gates) remain fixed and their connecting lines are redrawn, if needed.

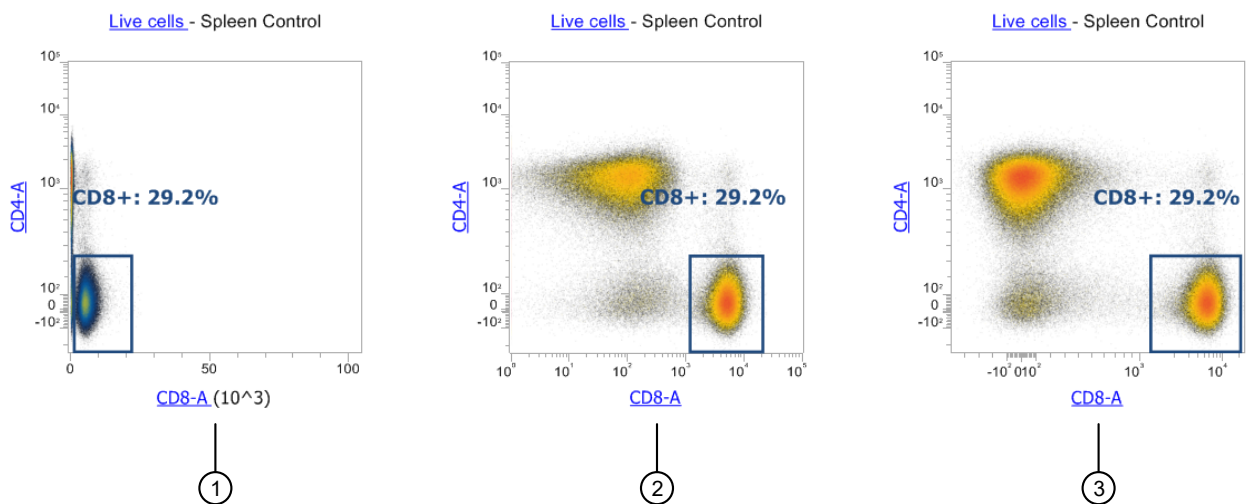


Figure 35 Changing plot scale type

- ① Linear
- ② Log
- ③ Hyperlog™

- For **Oval gates**, the coordinates of the bounding rectangle remain fixed and the perimeter ellipse and center position are redrawn, if needed.
- If the change in scale type results in a gate being off plot, the gate behaves as described in “Off-plot gates” on page 164.

## Changing plot parameter

- When the parameter of any axis is changed, including using the **Swap Axes** option in the **Plot context menu** (“Swap Axis” on page 192), all gates are maintained on the plot in their specified locations.
- If the change in parameter results in a change in scale type or scale range, the gates conform to the rules defined for changing the plot scale or scale range as described above.

## Changing plot type

- When changing from a single parameter to a dual parameter plot or vice versa, all gates are removed from the plots.
- When changing from any dual parameter plot another type of dual parameter plot, the gates are maintained at their specified locations.

# Workspace statistics

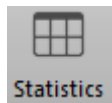
## Overview

The Attune™ Cytometric Software generates statistics from acquired events. Statistics can be displayed for any parameter and calculated for any defined population. As the events are acquired, statistics are updated in real time.

- To add a new **Statistics box** to the current **Workspace**, click the **Statistics** button on the **Workspace ribbon tab** (“Other group” on page 85).

Alternatively, select **Insert ▶ Statistics** from the

**Workspace view context menu** (“Workspace view context menu” on page 185) or the **Plot context menu** (“Plot context menu” on page 189).



Experiment: **In Vivo T cell Proliferation**  
Group: **Group**  
Sample: **Spleen Control**  
Time Recorded: **14:58:49**

Name	Gate	Count	%Total	%Gated
△ All Events	All Events	2,239,344	100.0	100.0
△ Cells	Cells	1,183,954	52.9	52.9
△ singlets	singlets	1,160,149	51.8	98.0
△ CD3+	CD3+	500,075	22.3	43.1
△ Live cells	Live cells	482,357	21.5	96.5
△ CD8+	CD8+	141,087	6.3	29.2
△ R7	R7	9	0.0	0.0
△ R8	R8	45	0.0	0.0

- The statistics available are: **Plate, Experiment, Group, Sample, Workspace, Plot Title, Gate, X parameter, Y parameter, Autogate Status, Comp Source, Count** (event count), **Events/μL, % Total, % Gated, Volume (μL), X Mean, Y Mean, X Median, Y Median, X Peak, Y Peak, X SD, Y SD, X %CV, Y %CV, X rSD, Y rSD, X %rCV, and Y %rCV.**
- You can customize the **Statistics box** using the **Customize Statistics** options available in the **Customize panel** (“Customize statistics box options” on page 445).
- When the statistics container is not large enough to hold all columns and rows of the **Statistics table**, indicators reveal that more data are present to the right or below the viewable area. To view any non-visible results, you must expand the results area.

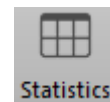
Experiment: **In Vivo T cell Proliferation**  
Group: **Group**  
Sample: **Spleen Control**  
Time Recorded: **14:58:49**

Name	Gate	Count	%Total	%Gated
△ All Events	All Events	2,239,344	100.0	100.0
△ Cells	Cells	1,183,954	52.9	52.9
△ singlets	singlets	1,160,149	51.8	98.0
△ CD3+	CD3+	500,075	22.3	43.1
△ Live cells	Live cells	482,357	21.5	96.5
△ CD8+	CD8+	141,087	6.3	29.2
△ R7	R7	9	0.0	0.0
△ R8	R8	45	0.0	0.0

- Double-clicking the **Statistics table** lets you scroll to see the contents that are outside the visible area. In this mode, the **right-click context menu** also includes an option to export the statistics of the selected table.

## Workspace statistics

- If no Workspace object is selected, clicking on the **Statistics** button on the **Workspace ribbon tab** or selecting **Insert > Statistics** from the **Workspace view context menu** inserts a **Workspace Statistics table** that shows the results from all gates.



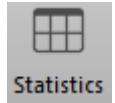
Experiment: **Whole Blood 1-4000 dilution**  
 Group: **Whole Blood**  
 Sample: **T6**  
 Time Recorded: **09:23:04**

Name	Count	Concentration	%Total	%Gated
△ All Events	57,545	30.4	100.0	100.0
△ PLT + WBC	57,545	30.4	100.0	100.0
Platelets	53,992	28.5	93.8	93.8
WBCs	3,430	1.8	6.0	6.0
-/+	84	0.0	0.1	0.1
+/+	2,456	1.3	4.3	4.3
-/-	53,893	28.5	93.7	93.7
+/-	1,112	0.6	1.9	1.9
RBC1	0	0.0	0.0	0.0
Noise	0	0.0	0.0	0.0

- The gates in the **Workspace Statistics table** are shown in the order specified by the **gate hierarchy**. Gates at the same hierarchical level are shown in the order they are created. The color of the gate is shown as a colored square to the left of the gate name.
- The **All Events** statistic node is shown at the top of the hierarchy. The **All Events** node in the **Statistics table** only shows results for these statistics: **%Gated**, **%Total**, and **Count**. All other statistics for the **All Events** node show **N/A**.

## Plot statistics (single plot)

- If a single plot is selected, clicking the **Statistics** button on the **Workspace ribbon tab** or selecting **Insert ▶ Statistics** from the **Plot context menu** creates a **Plot statistics box**, which displays the results from the gates present on the selected plot.



Parameters: **Blue 488 nm SSC - SSC-H vs Violet 405 nm SSC - VL1-H**  
Gate: **All Events**  
Experiment: **Whole Blood 1-4000 dilution**  
Group: **Whole Blood**  
Sample: **T6**  
Time Recorded: **09:23:04**

Name	Count	Concentration	%Total	%Gated
▾ All Events	57,545	30.4	100.0	100.0
▫ PLT + WBC	57,545	30.4	100.0	100.0
▫ RBC1	0	0.0	0.0	0.0
▫ Noise	0	0.0	0.0	0.0

- The **statistics box** for the selected plot also shows the **All Events** statistic node as well as the results from any gate present on the plot.
- The single **plot statistics box** shows the results for the gates on the current plot and any upstream gates.

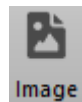
## Plot statistics (multiple plots)

- If multiple plots are selected, clicking the **Statistics** button on the **Workspace ribbon tab** or selecting **Insert ▶ Statistics** from the **Plot context menu** creates a **Plot statistic box** for each plot selected.
- Each **statistic box** contains the **All Events** statistics row, and only the statistics for the gates present on the selected plots as well as their parent gates.

## Workspace images and text

### Workspace images

You can add an **image** to the **Workspace** by selecting the **Image** button on the **Workspace ribbon tab** (“Other group” on page 85).



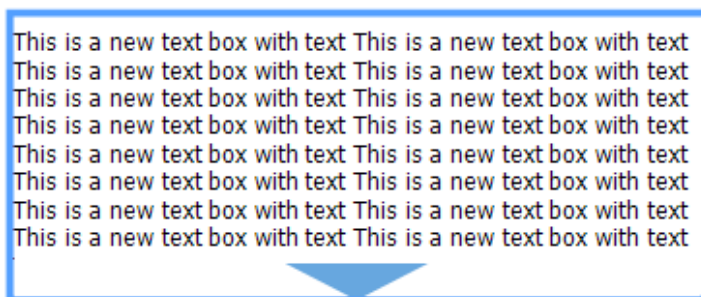
- The image types available for insertion include **JPG**, **GIF**, **BMP**, **PNG**, **TIF**, and **EMF** (Windows™ enhanced metafile).
- The image file cannot be larger than one page, and it is scaled to fit on the page and preserve the aspect ratio, if needed.
- Workspace images can be resized as described in “Resize Workspace objects” on page 123.

### Workspace text

You can add a **Text box** to the **Workspace** by selecting the **Text** button on the **Workspace ribbon tab** (“Other group” on page 85).



- The **Text box** displays the default text **New Text Box**, which can be edited in place without opening a separate dialog box. The **Text box** can contain at up to 500 characters.
- The style and border of the **Text box** can be customized as described in “Customize text box options” on page 444.
- If the **Text box container** is not large enough to hold all text, indicators reveal that more text is present to the right or below the viewable area. To view any non-visible text, you must expand the **Text box**.

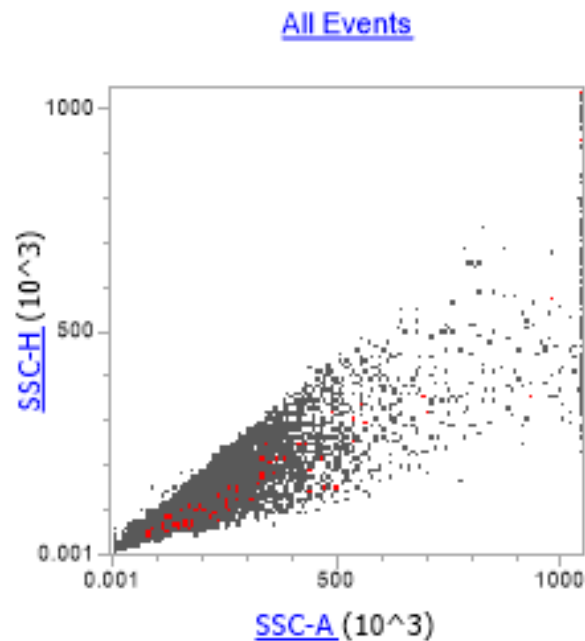




## Image backgating

**Image backgating** lets you identify the event that is associated with the captured cell image on a **Dot plot** or a **Precedence Density plot** in **Workspace view**.

When image backgating is applied, the events that have a corresponding image (**ImageFlag** parameter value of 1) are identified in the plot in red.



By default, backgated events appear on the associated plot in red, but you can select a different color for backgated events using the **Image Options dialog** (“Image Options” on page 673).

The backgated image events are always maintained as the top-most events in the precedence order of backgated events (i.e., they are painted on top of other events).

Image backgating persists at the Workspace level for an Experiment, Group, or Sample.

You can use the following methods to backgate images:

- Right-click the **cell image** you want to backgate in **Image View**, then select **Show Image on Plots**. You can select the Active image or any image from the **Image Gallery**.

To backgate multiple images, select the images you want to backgate in the **Image Gallery**, then right-click and select **Show Images on Plots**.

To undo the backgating of an image, right-click the backgated image, then select **Clear Image on Plots**.

To undo all backgates on a plot, right-click the Active image or on any image in the **Image Gallery**, then select **Clear All Images on Plots**.

- Click **Backgate All Images** in the **Image Settings ribbon tab** (“Image Settings tab” on page 112) to backgate all imaged events on the corresponding Workspace plots.

To undo the backgating of all images, click **Backgate All Images** again.

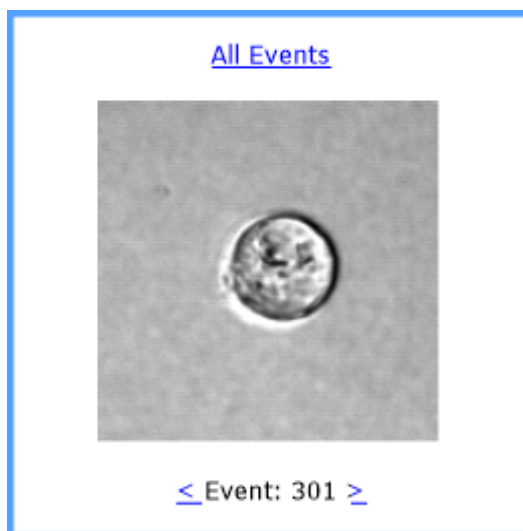
When enabled, the **Backgate All Images** button is shown in the **ON** state (blue). When disabled, the button is shown in its **OFF** state.

The **Backgate All Images** button is enabled only if the selected Sample’s listmode specification contains the **ImageFlag** and **Event** parameters.

## Cell image container

### Overview

**Cell Image Containers** enable you to add cell images specific to the sample to the current **Workspace**. They act like plots and can be batch printed with the **Workspace** (cell images cannot be printed from the **Image Gallery** in **Image view**).

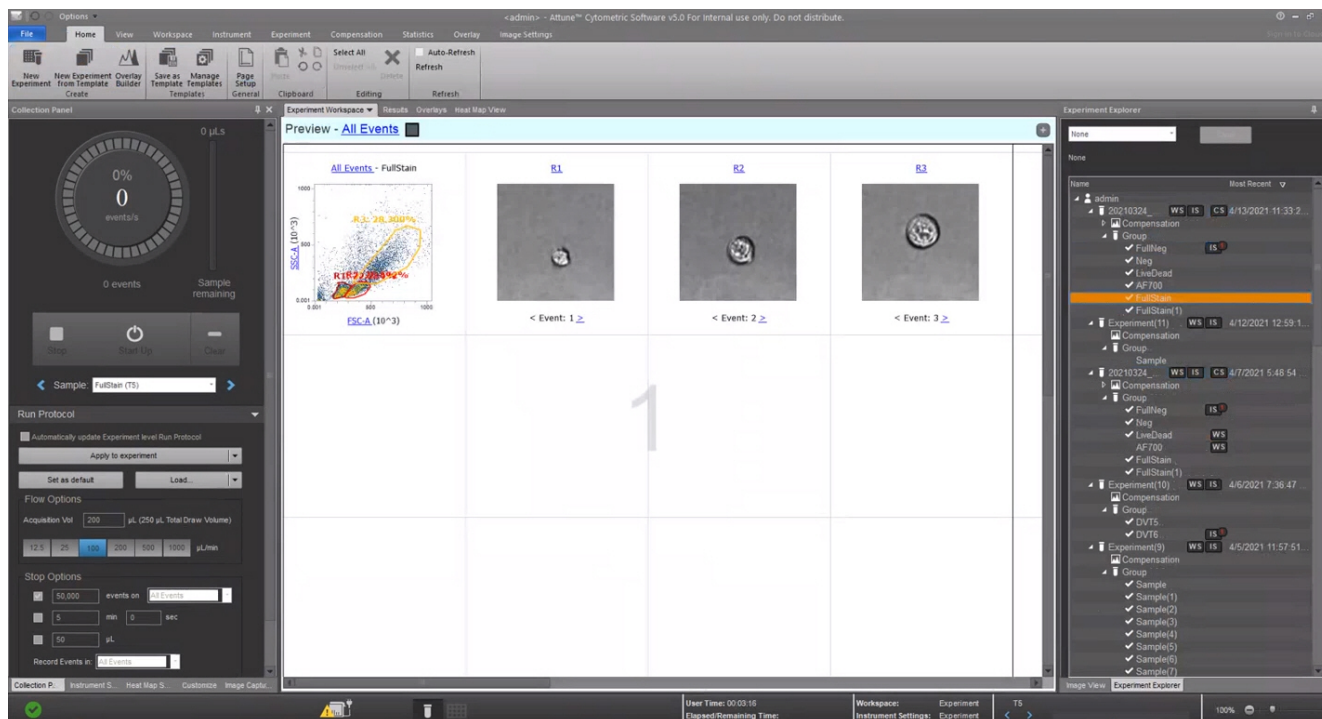


---

**Note:** **Cell image containers** are available only when using an Attune™ CytPix™ Flow Cytometer or if the active **Experiment** was created with an Attune™ CytPix™ Flow Cytometer.

---

In the following example, the **Workspace view** contains a **Dot plot** with three gated populations (**R1**, **R2**, **R3**) and three **cell image containers**. The **cell image containers** each have been gated to contain cell images from within one of the three gates (see “Population menu” on page 182).



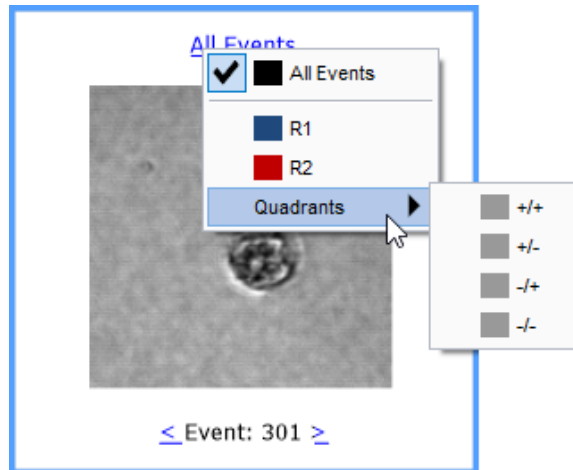
**Note:** For the **Cell Image Container context menu**, see “Cell Image Container context menu” on page 210.

## Population menu

- The **Cell Image Container** contains a **gate hyperlink** that lets you filter the images displayed in the container.  
By default, the gate hyperlink is set to **All Events**.



- Clicking the **gate hyperlink** opens the **Gate context menu** where you can select a gate to only allow images within the selected gate to be displayed.



- You can select any **gate type**. **Quadrant gates** are displayed in a submenu.
- The active gate is indicated by the check mark to the left of the gate name.
- When you select a gate, the cell image container only allows the selection of images where the events in the gate have a corresponding image.
- If the gate does not contain any images, the **cell image container** shows the message **No images are available**.
- When a gate is moved or the plot/gate is deleted, the **cell image container** updates accordingly. If the gate equation is changed, the **cell image container** can only show images within the gating hierarchy.

## Navigation

- The **Cell Image Container** shows the **event ID** as a label under the image.

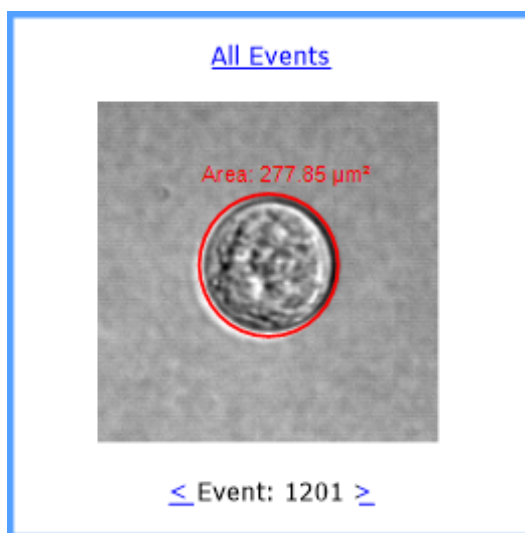


- The image in the **cell image container** can be changed in the selected gate using the < and > hyperlinks to the left and right of the event label that shows the active **event ID** of the image.

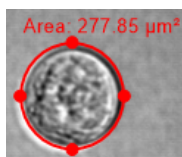
- The < hyperlink is disabled when the first available image in the selection is shown and the > hyperlink is disabled when the last image within the gate selection is displayed.
- The **selected images index** persists as part of the image container report object. The **image index** is the index of the image within a vector of images; it is **not** the **event ID**.
- If the **image index** does not exist when opening another sample, it defaults to the first image within the available images for the selected gate.

## Measure Image tool

The **Measure Image** tool enables you to measure the area of an imaged cell by drawing an ellipse in the **Cell Image Container** that encircles the imaged cell.



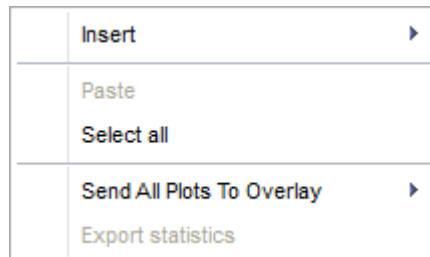
- When the **Measure** option is selected in the **Cell Image Container context menu** ("Cell Image Container context menu" on page 210) or the **Measure Image** tool is selected in **Cell Image Analysis** group of the **Workspace tab**, the measure image insert cursor is displayed.
- You can resize the ellipse using the handles that become visible when the ellipse is selected to stretch and rotate it.
- To delete the measurement tool, click the ellipse, then select from the keyboard or the context menu.
- The measured area is based on a conversion factor of micrometers per pixel.



## Workspace view context menu

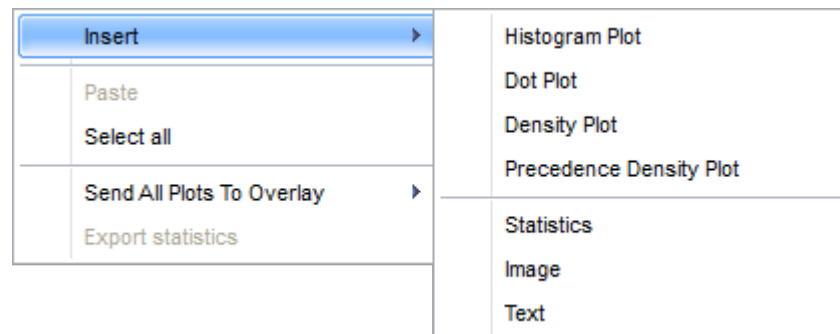
### Overview

The **Workspace view context menu** is displayed when you right-click any empty area (i.e., white space) within the **Workspace view**. It contains the tools to insert plots, statistics, images, and text to the Workspace, to paste and select Workspace objects, and to export statistics.



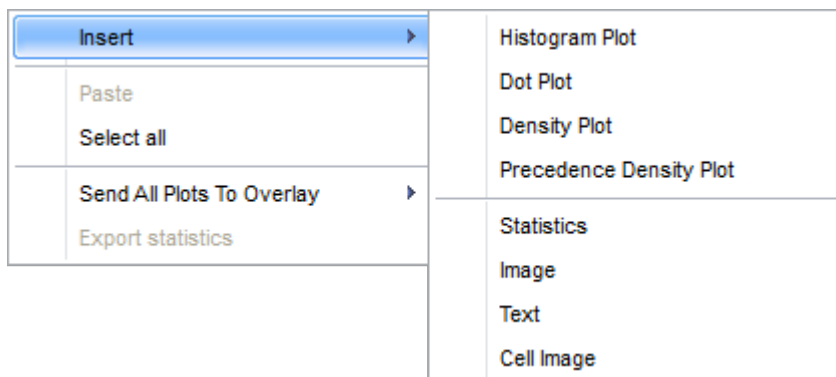
### Insert

Selecting the **Insert** option displays a submenu, allowing you to add **Histogram Plot**, **Dot Plot**, **Density Plot**, **Precedence Density Plot**, **Statistics**, **Image**, and **Text** objects to the current **Workspace**.



These menu options have the same effect as pressing the respective buttons on the **Workspace ribbon tab** as described on “Workspace tab” on page 81.

If you are using an Attune™ CytPix™ Flow Cytometer or when an **Experiment** that was created with an Attune™ CytPix™ Flow Cytometer is active, the Insert submenu also includes an option to insert **Cell Image Container** to the current **Workspace**.



## Paste and Select all

### Paste

- **Paste** inserts the object currently in the clipboard to the Workspace at the location clicked. Paste option is only available if a Workspace object has been copied or cut to the clipboard.
- Clicking the **Ctrl+V** keyboard combination pastes the object at same location as the original. As each new object is pasted, it is inserted cascading downwards and to the right of the original location.
- Objects can be cut or copied from any Workspace and pasted to any other Workspace.
- Plots are pasted with any gates present. If a gated plot is copied to a new location where the parent gate does not exist, the pasted plot will not be gated and will display all events.
- If a Plot statistics box is pasted to a new Workspace and the target Workspace does not contain a copy of the original plot, the statistics box will update to show all gates present in the current Workspace. The Plot Axis and Gate fields will be removed from the statistics box header (if originally present).
- Pasted objects are selected by default after the paste action has completed.

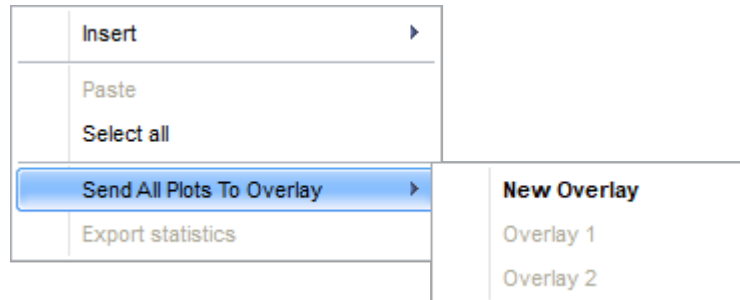


## Select all

- **Select all** selects all objects on the Workspace.

## Send All Plots To Overlay

- **Send All Plots to Overlay** sends all Workspace plots to a selected **Overlay** (Chapter 9, “Overlays”).

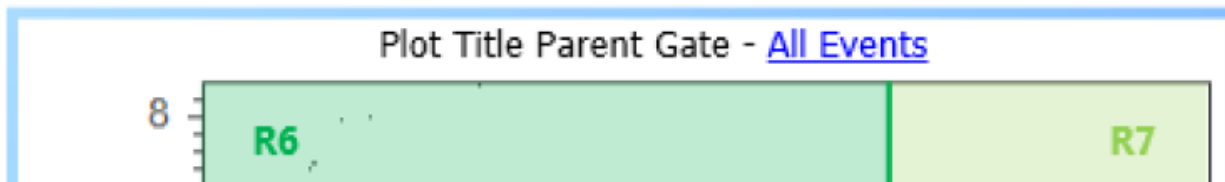


The options depend on the type of plot present on the **Workspace** and the **Overlay plots** already present in the **Overlay view**.

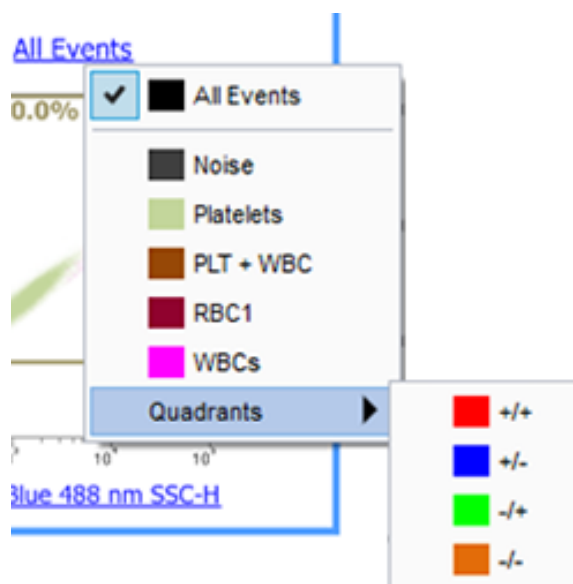
- The **New Overlay** option is always present and enabled unless the maximum number of **Overlays** have already been created.
- The names of any **Overlay plots** already present in the **Overlay** application are also displayed.
- If both **Histogram** and dual parameter plots (**Dot plot** or **Density plot**) are present on the **Workspace**, the only option is to send plots to a **New Overlay**. This action sends Histogram plots to a new Overlay and dual parameter plots to another new Overlay.
- If only **Histogram plots** or dual parameter plots (**Dot plot** or **Density plot**) are present on the **Workspace**, you can either create a **New Overlay** or send the plots to an existing **Overlay** that contains plots with the same number of parameters.  
**Overlay plot** names in the context menu are disabled for **Overlay plots** with non-matching number of parameters.
- The **Overlay name** in the context menu is also disabled if the number of plots to be sent exceeds the allowable number of plots in the specific **Overlay**.

## Plot title hyperlink

If **Include** option is selected in the **Customize Plot** options (“Customize plot options” on page 421), the plot title displays the current parent gate of the plot as a **hyperlink**. This option is selected by default.



- Clicking the **plot title hyperlink** displays a dropdown menu that contains a list of all available gates, including any derived gates.  
Any gates present on the currently selected plot and their daughter gates are not included in the list of gates.  
Quadrant gates are displayed in a submenu.



- You can select one gate from the dropdown menu to be used as the parent gate of the current plot. The hyperlink then updates to show the currently selected gate name.

# Plot context menu

## Overview

The **Plot context menu** is opened when you right-click any empty area (i.e., white space) within the plot boundary, but not on another active area (gate, plot title, or plot axis).

The **Plot context menu** contains the tools for gate creation and customization, to cut, copy, paste, and delete selected items, and save all selected items as a new graphics file.

Customize
Scale ▶
Set Population ▶
Add gate ▶
Insert plot statistics
Swap Axis
Cut
Copy
Paste
Send to overlay ▶
Delete
Save as

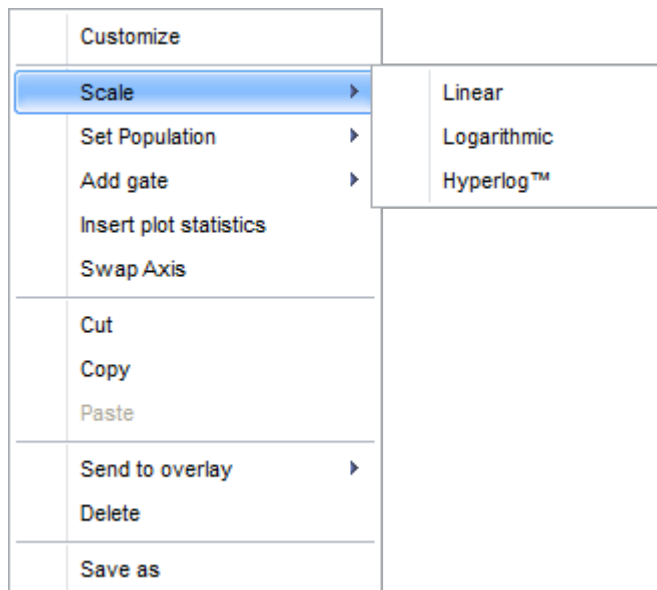
## Customize

Selecting **Customize** from the dropdown menu displays the **Plot Customize panel**, which enables you to format different properties of the selected plot as described in “Customize plot options” on page 421.

- If the panel is closed, it will be opened and positioned in the left docking panel.
- If the panel is docked and not in the foreground, it will be brought to the front of the docked panels.

## Scale

**Scale** sets all parameter axes to the selected scale. Available scale types are **Linear**, **Logarithmic**, and **Hyperlog™**.



- The **Scale** option is available only when single or multiple plots are selected. If images or statistics objects are also selected, this option is not available.
- On dual-parameter plots (**Dot** or **Density plots**) the X and Y-axes are modified.  
On single parameter plots (i.e., **Histograms**), the selection is only applied to the X-axis.

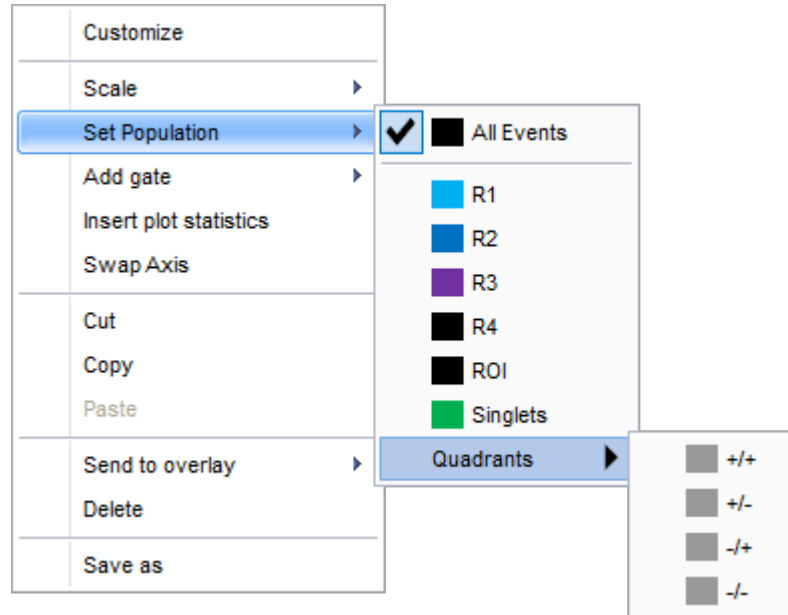
---

**Note:** The **Hyperlog™** scale uses log-linear hybrid transformations to show compensated flow cytometry data that frequently contain negative values due to compensation. Logarithmic transformations cannot properly handle negative values, and poorly show normally distributed cell types.

---

## Set Population

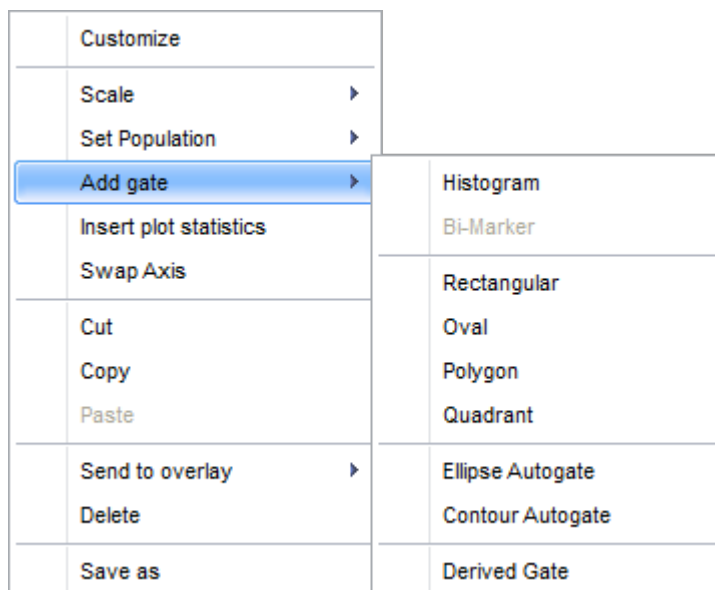
**Set Population** lets you limit the data shown on a plot to a given gate. It makes the plot a **daughter plot** of the selected **upstream gate**.



- A single gate can be selected from the list of available gates to be used as a **parent gate** for the selected plot. The selected gate is indicated by a tick mark in the list. It makes the plot a **daughter plot** of the **upstream gate**.
- The list of gates includes the **All Events** option and any available derived gates.
- The list does not include any gates present on the selected plot or any gates derived from those gates.
- The **Set Population** option is also available when multiple plots are selected. In such cases, the available gate list excludes the gates that are present on any of the selected plots or any gates derived from those gates. The selected gate is used as a parent gate for all selected plots.

## Add gate

**Add gate** enables you to create a new gate or a new derived gate in the default position on the currently selected plot. This option is available only when a single plot is selected.



- The **Histogram** option is only available for **Histogram plots**.  
**Rectangular**, **Oval**, **Polygon**, **Ellipse Autogate**, **Contour Autogate**, and **Quadrant** options are available only for dual parameter plots.  
Selecting these options creates the selected gate type in the default position described in “Create and delete gates” on page 145.
- The **Derived Gate** option opens the **Derived Gate dialog** (see “Derived gate dialog” on page 739).

## Insert plot statistics

**Insert plot statistics** inserts a **Plot statistics box** for each selected plot as described on “Plot statistics (single plot)” on page 177. This option is available when single or multiple plots are selected.

## Swap Axis

**Swap Axis** flips the axes of the selected plot, so that the X parameter becomes the Y parameter, and the Y parameter becomes the X parameter.

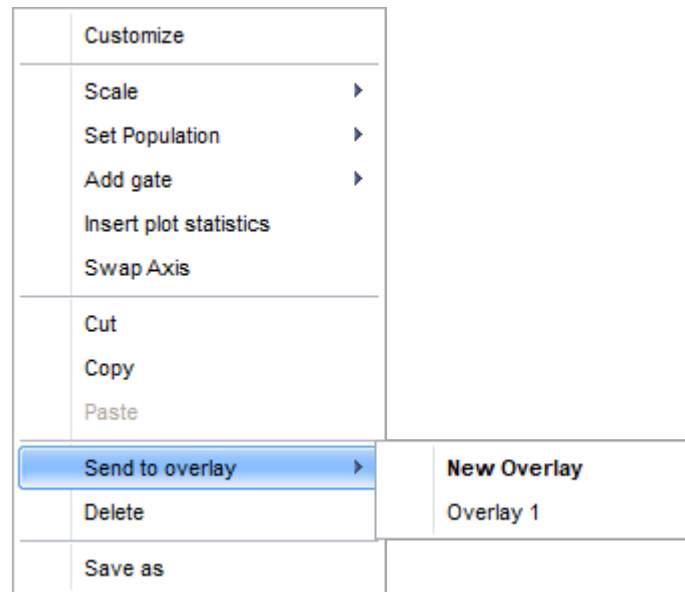
The **Swap Axis** option is only visible if at least one dual parameter plot is selected.

## Cut, Copy, and Delete

- **Cut** copies the selected items to the clipboard and removes them from the **Workspace**. Cut items can be pasted as described in “Paste Workspace objects” on page 125.
- **Copy** copies the selected items to the clipboard, but keeps the original items on the **Workspace**. Copied items can be pasted as described in “Paste Workspace objects” on page 125.
- **Delete** permanently removes all selected items from the **Workspace**.

## Send to overlay

**Send to overlay** sends all selected plots to a selected **Overlay** (Chapter 9, “Overlays”). The available options depend on the type of plots selected and the **Overlay plots** already present in the **Overlay view**.



- The **New Overlay** option is always present and enabled unless the maximum number of **Overlays** have already been created.
- The names of any **Overlay plots** already present in the **Overlay application** are also displayed.
- If both **Histogram** and dual parameter plots (**Dot plot** or **Density plot**) are selected, the only option is to send plots to a **New Overlay**. This action sends histogram plots to a new Overlay and dual parameter plots to another new Overlay.
- If only **Histogram** plots are present or only dual parameter plots are selected, you can either create a **New Overlay** or send the plots to an **existing Overlay** that contains plots with the same number of parameters.
- **Overlay plot names** in the context menu are disabled for **Overlay plots** with non-matching number of parameters. The **Overlay name** in the context menu is also disabled if the number of plots to be sent exceeds the allowable number of plots in the specific Overlay.

## Save as

**Save as** opens the **File Save (Export) dialog** as described in “File Save (Export) dialog” on page 715, which lets you send all selected items to a new graphics file.

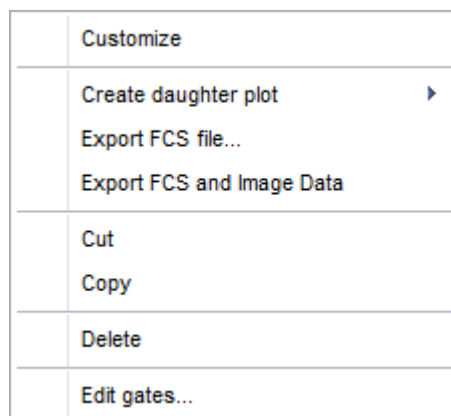
- The layout of the graphics file replicates the layout of the selected items on the current **Workspace**.
- The selection area for saving can only cover one page.
- The available image file formats are **PDF, BMP, JPG, TIFF, GIF, PNG**, and **EMF** (Windows™ Metafile). The default format is **PNG**.

## Gate context menu

### Overview

The **Gate context menu** is displayed when you right-click any part of a **Gate**, the **gate boundary line**, or a **control point**.

The **Gate context menu** contains the tools for customizing and editing gates, creating daughter plots, exporting FCS file, exporting FCS and image data, and for cutting, copying, deleting, and editing selected gates.



### Customize

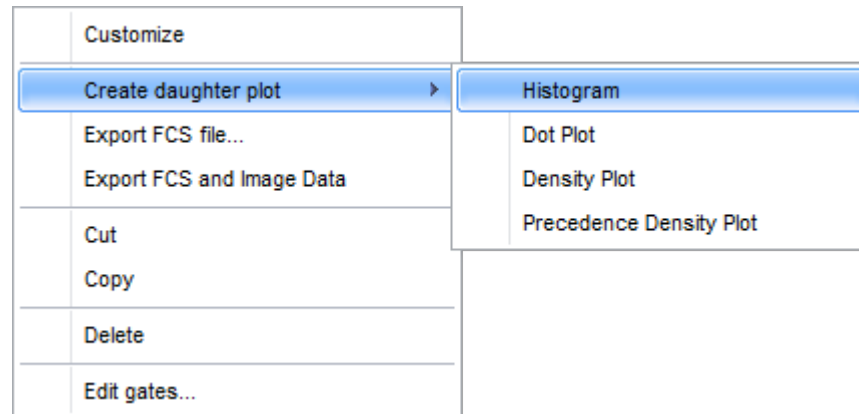
**Customize** shows the **Gate Customize panel**, which enables you to format different properties of the selected gate as described in “Customize gate options” on page 447.

- If the panel is closed, it will be opened and positioned in the left docking panel.
- If the panel is docked and not in the foreground, it will be brought to the front of the docked panels.



## Create daughter plot

**Create daughter plot** opens a submenu, which enables you to select the type of plot to create. This creates a new plot gated on the currently selected gate.



- Properties of the newly created **daughter plot** are based on the defaults set in the **Plot Options dialog** (“Fonts and Styles” on page 649) and not those of the **parent plot**.
- If you select a dual parameter plot type (**Dot plot**, **Density plot**, or **Precedence Density plot**), the new plot displays parameter 1 on the X-axis (normally FSC) and parameter 2 (normally SSC) on the Y-axis. If FSC and/or SSC are not being collected, the appropriate parameter 1 and 2 are used as described in “Plots” on page 138.
- If you select the **Histogram** option, the new plot displays parameter 1 (normally FSC) on the X-axis and Count on the Y-axis.
- If the **Workspace** is in **Freeform** mode (“Freeform mode” on page 126), the new plot is created in the default position as described ““Create Workspace objects” on page 120”.
- If the **Workspace** is in **Auto Layout** mode (“Auto Layout mode” on page 127), a new plot of the same size as the parent is inserted in the next free grid location available on the **Workspace**.

## Export FCS file

**Export FCS file** opens the **File Save dialog** to save a new FCS file containing only the events within the selected gate.

- This option is only enabled if the active sample has data and acquisition is not in progress.
- This option is inactive for quadrants.
- If there are any values for any of the keywords that have changed, a dialog is displayed allowing you to update the keywords.
  - If you select **Ignore**, the keywords are not updated (except for originality keywords).
  - If you select **Cancel**, the gated FCS data are not exported.
  - If you select **Update keywords** option, the keywords that have changed are updated.

- In all cases, the following keywords are updated:
  - **\$ORIGINALITY**: Set to NonDataModified
  - **\$LAST\_MODIFIED**: The time a data file is appended or updated.
  - **\$LAST\_MODIFIER**: The name of the currently logged in user.
  - **\$TOT**: Updated to reflect the number of events exported in the gate.
- When exporting an FCS file, the exported data can be appended to the original FCS file, which creates a single FCS file that includes the new data from the gate.  
To add the exported data can be appended to the original FCS file, select **Extend FCS File** on the **File Save dialog**. This option can be used for whole datasets or specifically for exporting from a gate.  
When **Extend FCS File** option is selected, the original FCS file is appended with the new data, and the **\$ORIGINALITY**, **\$LAST\_MODIFIER**, and **\$LAST\_MODIFIED** keywords in the FCS file are updated to indicate that the data have been appended and modified.

## Cut, Copy, and Delete

### Cut

**Cut** copies the selected gate to the clipboard and removes it from the plot.

When a gate is removed from a plot, any Derived Gate equations that include the removed gate are modified to exclude that gate. If the Derived gate includes only the gate that was removed, then the entire Derived Gate equation is deleted.

The cut gate can be pasted as described on “Paste Workspace objects” on page 125.

### Copy

**Copy** copies the selected gate to the clipboard, but keeps the original gate on the plot.

### Delete

**Delete** permanently deletes the currently selected gate and removes the gate from any Derived Gate equations and plot gates. If the gate equation includes only the deleted gate, then the entire gate equation is deleted.

## Edit gates

**Edit gates** opens the **Edit Gates dialog** (“Edit gates dialog” on page 742).

The **Edit Gates dialog** provides a list of all available gates on the active **Workspace** and enables you to edit the gate color, gate math expression, and the order in which the gates are displayed in the dialog, to backgate all plots, and to delete gates.

# Plot axis context menu

## Overview

The **Plot axis context menu** is displayed when you right click a **plot scale** or a **parameter name**.

The **Plot axis context menu** contains the tools for creating and customizing plots and gates, to cut, copy, paste, and delete selected items, and save all selected items as a new graphics file.

Customize
Scale ▶
Set Population ▶
Backgates ▶
Add gate ▶
Insert plot statistics
Swap Axis
Cut
Copy
Paste
Send to overlay ▶
Delete
Save as

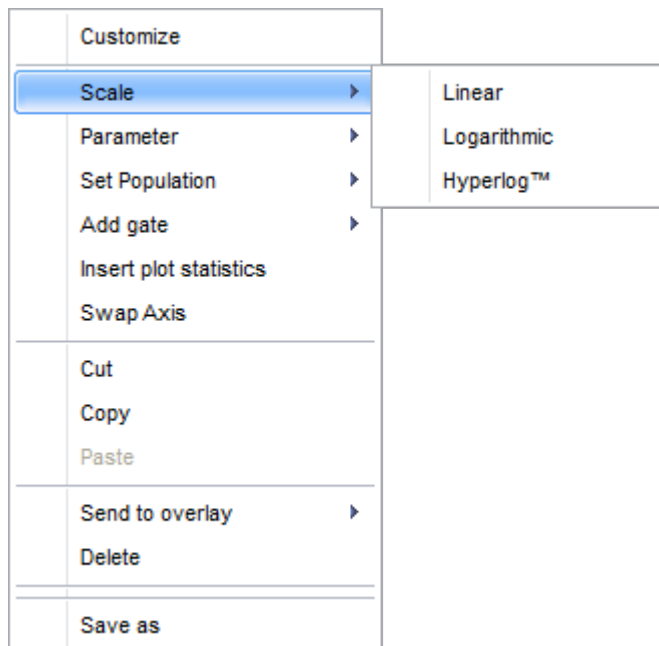
## Customize

**Customize** displays the **Plot Customize panel**, which enables you to format different properties of the selected plot as described in “Customize plot options” on page 421.

- If the panel is closed, it will be opened and positioned in the left docking panel.
- If the panel is docked and not in the foreground, it will be brought to the front of the docked panels.

## Scale

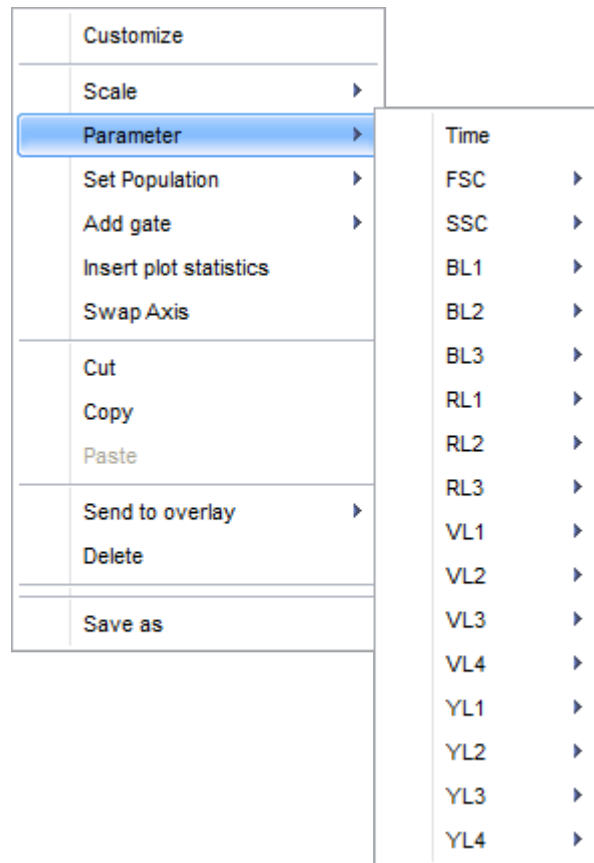
**Scale** sets all parameter axes to the selected scale. Available scale types are **Linear**, **Logarithmic**, and **Hyperlog™**.



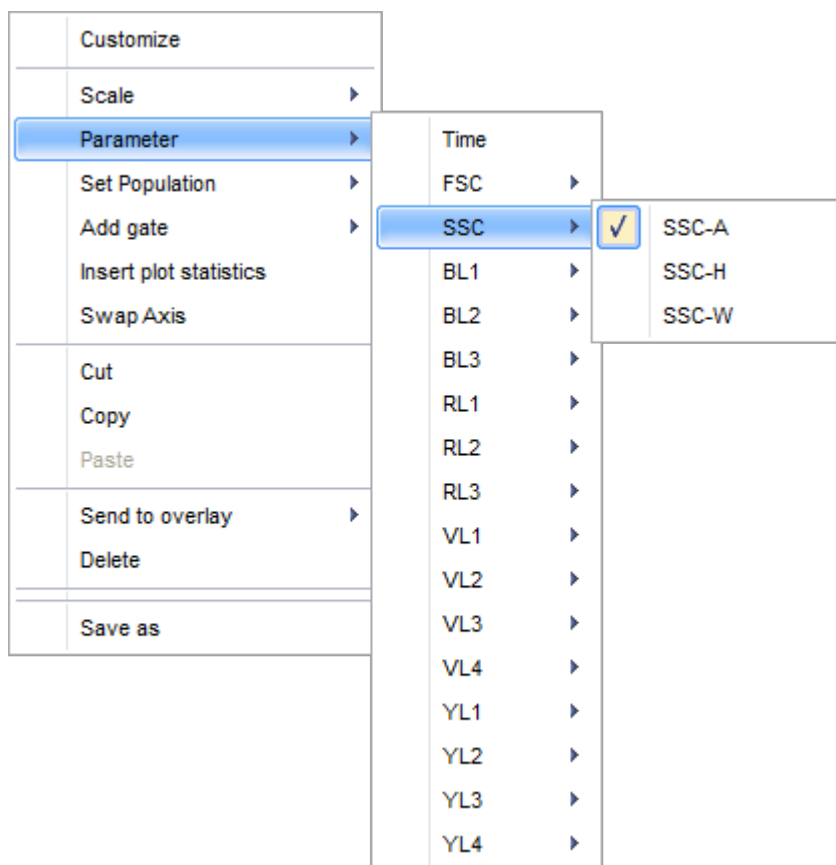
- The **Scale** option is available only when single or multiple plots are selected. If images or statistics objects are also selected, this option is not available.
- On dual-parameter plots (**Dot plots** or **Density plots**) the X- and Y-axes are modified. On single parameter plots (i.e., **Histograms**), the selection is only applied to the X-axis.

## Parameter

**Parameter** lets you select the parameters to show on a selected plot axis. This option is visible and enabled when single or multiple plots are selected. If images or statistics objects are also selected, this option is not visible.

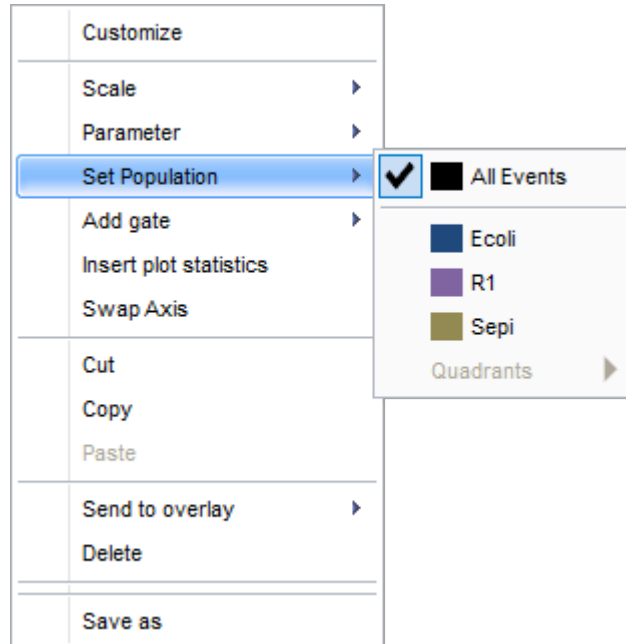


- The options displayed in the **Parameter submenu** depend on the number of parameters present in the current sample.  
If no dataset is present (i.e., the sample is empty), then the parameter list is based on the available and enabled parameters in the **Instrument Settings** (“Parameters” on page 385).
- If 18 or less parameters are selected in the current sample, the **Parameter submenu** displays all the parameters.
- The currently shown parameter appears checked in the menu, and you can select any of the available parameters to show. The appropriate plot axis is updated on all selected plots to show the chosen parameter.
- If more than 18 parameters are selected in the current sample, the **Plot axis context menu** displays the parameter names with the parameter type displayed as a submenu.



## Set Population

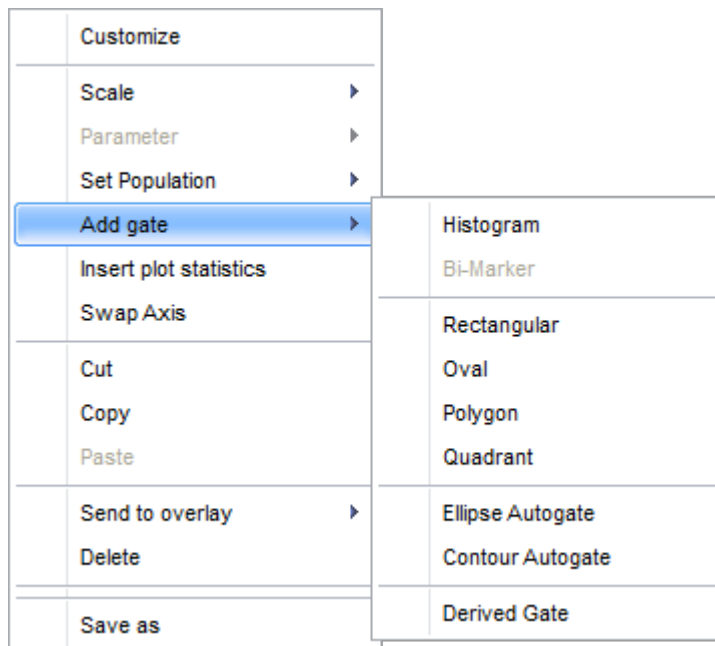
**Set Population** enables you to limit the data displayed on a plot to a given gate. It makes the plot a **daughter plot** of the selected **upstream gate**.



- A single gate can be selected from the list of available gates to be used as a **parent gate** for the selected plot. The selected gate is indicated by a tick mark in the list. It makes the plot a **daughter** of the **upstream gate**.
- The list of gates includes the **All Events** option and any available derived gates.
- The list does not include any gates present on the selected plot or any gates derived from those gates.
- The **Set Population** option is also available when multiple plots are selected. In such cases, the available gate list excludes the gates that are present on any of the selected plots or any gates derived from those gates. The selected gate is used as a parent gate for all selected plots.

## Add gate

**Add gate** enables you to create a new gate or a new derived gate in the default position on the currently selected plot. This option is available only when a single plot is selected.



- The **Histogram** option is only available for **Histogram plots**. **Rectangular**, **Oval**, **Polygon**, **Ellipse Autogate**, **Contour Autogate**, and **Quadrant** options are available only for dual parameter plots. Selecting these options creates the selected gate type in the default position described in “Create and delete gates” on page 145.
- The **Derived Gate** option opens the **Derived Gate dialog** described on “Derived gate dialog” on page 739.

## Insert plot statistics

**Insert plot statistics** inserts a **Plot statistics box** for each selected plot as described in “Plot statistics (single plot)” on page 177. This option is available when single or multiple plots are selected.

## Swap Axis

**Swap Axis** flips the axes of the selected plot, so that the X parameter becomes the Y parameter, and the Y parameter becomes the X parameter.

The **Swap Axis** option is only visible if at least one dual parameter plot is selected.

## Cut, Copy, Paste, and Delete

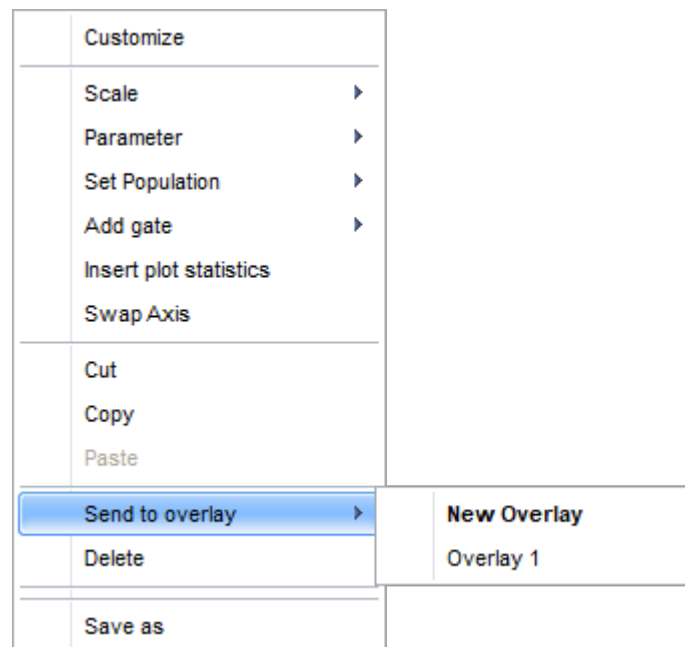
- **Cut** copies the selected items to the clipboard and removes them from the **Workspace**.
- **Copy** copies the selected items to the clipboard, but keeps the original items on the **Workspace**.



- **Paste** inserts the gate currently in the clipboard to the **Workspace**. The new gate is assigned the next available gate name as described in “Gate name” on page 146.
  - The size and position of the pasted gate will be identical to the location and size of the original gate when it was copied to the clipboard, even if the location is outside the current plot area.
  - The **Paste** option is available only if a single plot is selected and if a gate type relevant to the current plot type has previously been copied into the clipboard.
  - If **Smart gating** is enabled, the new gate name is created as described on “Name options” on page 459.
- **Delete** permanently removes all selected items from the **Workspace**.

## Send to overlay

**Send to overlay** sends all selected plots to a selected **Overlay** (page 234). The available options depend on the type of plots selected and the **Overlay plots** already present in the **Overlay view**.



- The **New Overlay** option is always present and enabled unless the maximum number of **Overlays** have already been created.
- The names of any **Overlay plots** already present in the **Overlay application** are also displayed.
- If both **Histogram** and dual parameter plots (**Dot plot** or **Density plot**) are selected, the only option is to send plots to a **New Overlay**. This action sends histogram plots to a new Overlay and dual parameter plots to another new Overlay.
- If only **Histogram** plots are present or only dual parameter plots are selected, you can either create a **New Overlay** or send the plots to an **existing Overlay** that contains plots with the same number of parameters.
- **Overlay plot names** in the context menu are disabled for **Overlay plots** with non-matching number of parameters. The **Overlay name** in the context menu is also disabled if the number of plots to be sent exceeds the allowable number of plots in the specific Overlay.

## Save as

**Save as** opens the **File Save (Export) dialog** as described in “File Save (Export) dialog” on page 715, which lets you send all selected items to a new graphics file.

- The layout of the graphics file replicates the layout of the selected items on the current **Workspace**.
- The selection area for saving can only cover one page.
- The available image file formats are **PDF**, **BMP**, **JPG**, **TIFF**, **GIF**, **PNG**, and **EMF** (Windows™ Metafile). The default format is **PNG**.

## Plot parameter hyperlink

The **plot parameter name**, the **X-axis** and **Y-axis** of a dual parameter plot, or the **X-axis** of a **Histogram plot** are **hyperlinks**.

- Clicking on a **parameter name** displays the list of available parameters and calculated parameters. You select a parameter to apply to the selected axis from this list.
- The options shown in the **parameter name context menu** depend on the number of parameters present in the current sample. Only parameters that are part of the dataset and selected in the **Workspace filters** (“Workspace filter dropdown” on page 105) are listed.  
If no dataset is present (that is, the sample is empty), then the parameter list is based on the available parameters in the **Instrument Settings** (“Parameters” on page 385).
- If 18 or fewer parameters are selected in the current sample, the **Parameter submenu** shows all the parameters present in the current sample.

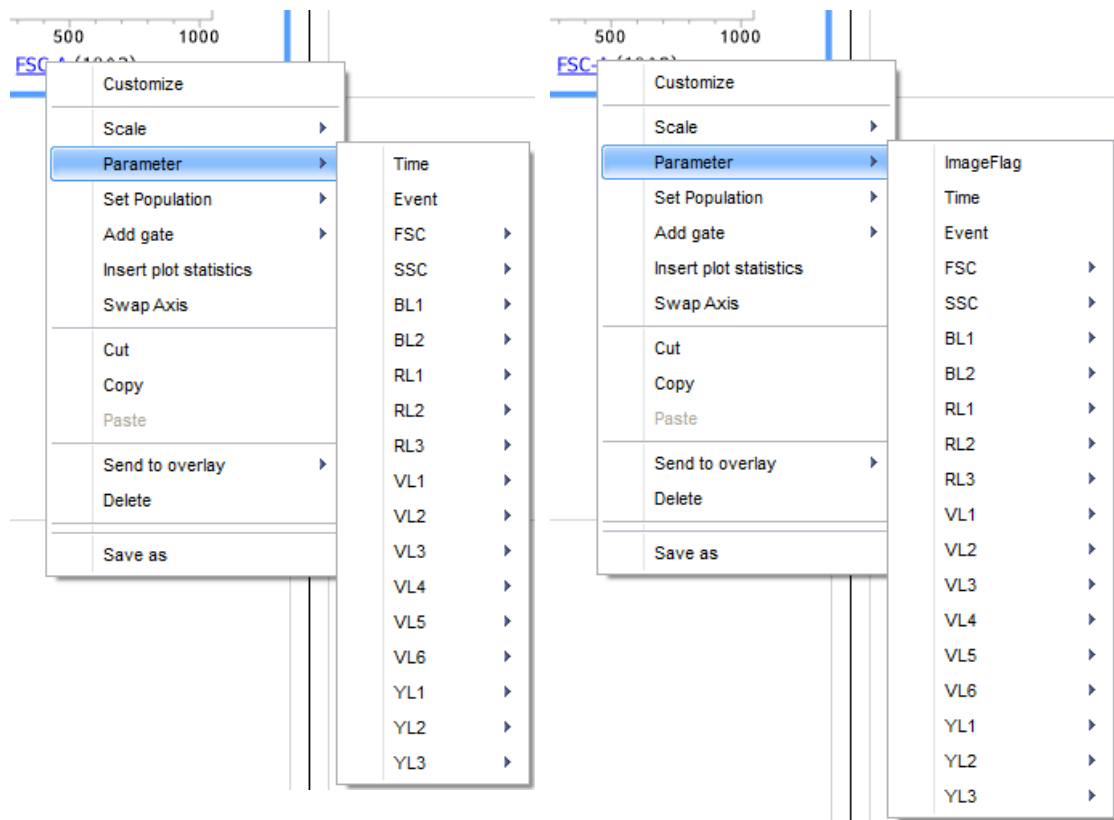
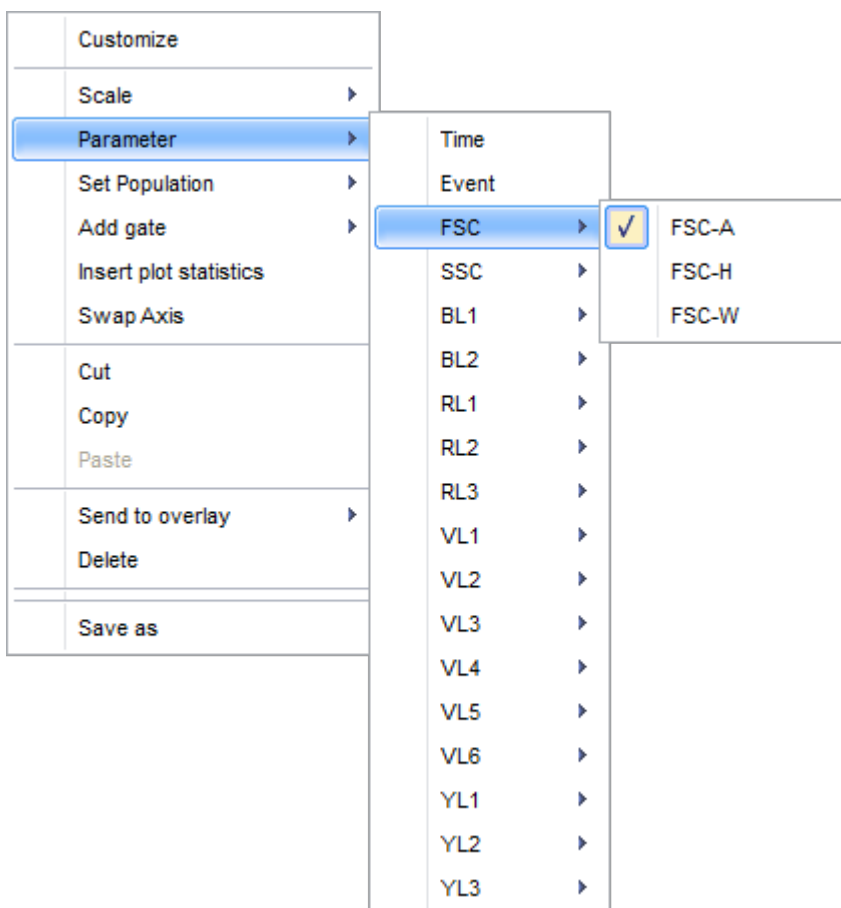


Figure 36 Parameter submenu in Attune™ NxT (left) and Attune™ CytPix™ Flow Cytometer(right). The list for the Attune™ CytPix™ Flow Cytometer shows the ImageFlag parameter only available in that instrument.

- The currently shown parameter appears checked in the menu, and you can select any of the available parameters to show. The appropriate plot axis is updated on all selected plots to show the selected parameter.



- If more than 18 parameters are selected in the current sample, the parameter name context menu displays the parameter names with the parameter type displayed as a submenu.
- If only one parameter type exists, the parameter choice remains in the Main menu.
- If multiple plots are selected, the chosen parameter is applied to the selected axis of all selected plots. Any Y-parameter choice is not applied to any Histogram plots that are part of the multiselect.

# Statistics context menu

## Overview

The **Statistics context menu** is shown when you right-click a **Statistics box** on the **Workspace** (“Workspace statistics” on page 175).

The **Statistics context menu** contains the tools for customizing the **Statistics box**, to cut, copy, and delete the selected **Statistics box**, and to export the selected **Statistics** as a single file.

	Customize
	Cut
	Copy
	Delete
	Export stats

## Customize

**Customize** displays the *Statistics Box Customize panel*, which allows you to customize the statistic style and formatting of the selected Statistics box as described in “Customize Statistics Box Options” on “Customize statistics box options” on page 445.

- If the panel is closed, it will be opened and positioned in the left docking panel.
- If the panel is docked and not in the foreground, it will be brought to the front of the docked panels.

## Cut, Copy, and Delete

- **Cut** copies the selected items to the clipboard and removes them from the **Workspace**. Cut items can be pasted as described in “Paste Workspace objects” on page 125.
- **Copy** copies the selected items to the clipboard, but keeps the original items on the **Workspace**. Copied items can be pasted as described in “Paste Workspace objects” on page 125.
- **Delete** permanently removes all selected items from the **Workspace**.

## Export stats

**Export stats** opens the **Save As dialog**, which lets you specify a file name and location to export the selected statistics table as a CSV file.

This option is only enabled if the statistics table is activated by double clicking the table.

## Image context menu

### Overview

The **Image context menu** is displayed when you right-click a **Workspace image** on the **Workspace** (“Workspace images” on page 178).

The **Image context menu** contains the tools to cut, copy, and delete the selected image.

	Cut
	Copy
	Delete

### Cut, Copy, and Delete

- **Cut** copies the selected images to the clipboard and removes them from the **Workspace**. Cut images can be pasted as described in “Paste Workspace objects” on page 125.
- **Copy** copies the selected images to the clipboard, but keeps the original images on the **Workspace**. Copied images can be pasted as described in “Paste Workspace objects” on page 125.
- **Delete** permanently removes all selected images from the **Workspace**.

## Text box context menu

### Overview

The **Text box context menu** is displayed when you right-click a **Text box** on the **Workspace** (“Workspace images and text” on page 178).

The **Text box context menu** contains the tools to customize, cut, copy, and delete the selected **Text box**.

	Customize
	Cut
	Copy
	Delete

## Customize

**Customize** displays the **Customize Text Box panel**, which enables you to change the font and style options of the text and border width of the selected **Text Box** as described in “Customize text box options” on page 444.

- If the panel is closed, it will be opened and positioned in the left docking panel.
- If the panel is docked and not in the foreground, it will be brought to the front of the docked panels.

## Cut, Copy, and Delete

- **Cut** copies the selected **Text box** to the clipboard and removes it from the **Workspace**. Cut **Text box** can be pasted as described in “Paste Workspace objects” on page 125.
- **Copy** copies the selected **Text box** to the clipboard, but keeps the original **Text box** on the **Workspace**. Copied **Text box** can be pasted as described in “Paste Workspace objects” on page 125.
- **Delete** permanently removes all selected **Text boxes** from the **Workspace**.

## Text box text context menu

### Overview

The **Text box text context menu** is displayed when you right-click the **text** within the edit area of a **Text box** while the **Text box** is being edited or the **text** selected.

The **Text box text context menu** contains the tools to cut, copy, and delete the selected text.

Cut	Ctrl+X
Copy	Ctrl+C
Paste	Ctrl+V

## Cut, Copy, and Delete

- **Cut** copies the selected text to the clipboard and removes it from the **Text box**. Cut text can be pasted as described in “Paste Workspace objects” on page 125.
- **Copy** copies the selected text to the clipboard, but keeps the original text on the **Text box**. Copied text can be pasted as described in “Paste Workspace objects” on page 125.
- **Delete** permanently removes all selected text from the **Text box**.

## Cell Image Container context menu

### Overview

The **Cell Image Container context menu** is displayed when you right-click a **Cell Image Container**.

The **Cell Image Container context menu** contains the tools to customize, cut, copy, delete, and measure the selected cell image.

Customize
Cut
Copy
Delete
Measure

### Customize

**Customize** displays the **Customize Image panel**, which enables you to change **Image Adjustment** and **Mask Settings** of the image in the selected Cell Image Container as described in “Customize Image options” on page 484.

- If the panel is closed, it will be opened and positioned in the left docking panel.
- If the panel is docked and not in the foreground, it will be brought to the front of the docked panels.

### Cut, Copy, and Delete

- **Cut** copies the selected **Cell Image Container** to the clipboard and removes it from the **Workspace**. Cut **Cell Image Container** can be pasted as described in “Paste Workspace objects” on page 125.
- **Copy** copies the selected **Cell Image Container** to the clipboard, but keeps the original **Cell Image Container** on the **Workspace**. Copied **Cell Image Container** can be pasted as described in “Paste Workspace objects” on page 125.
- **Delete** permanently removes all selected **Cell Image Containers** from the **Workspace**.

### Measure

**Measure** lets you draw a **measurement ellipse** in a **Cell Image Container** to measure the area of the cell in the image (“Measure Image tool” on page 184).

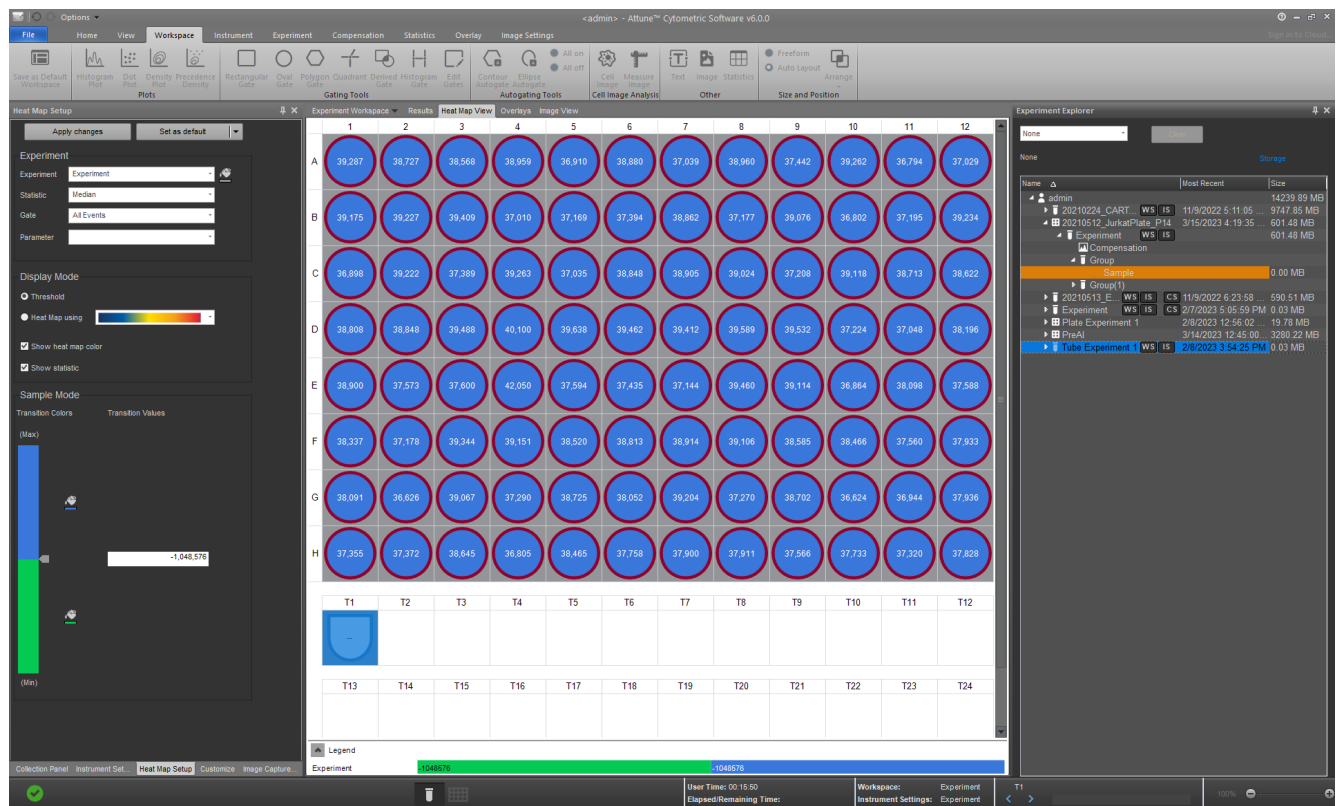




# Heat Map View

## Overview

The **Heat Map View** provides a graphical method for setting up and analyzing plate- and tube-based experiments. By default, the **Heat Map View** is docked to the **Main Application Workspace** (“Main application workspace” on page 56).



## Display of samples

### Overview

- The **Heat Map** displays both **Well** and **Tube Samples**.
- A tube- or plate-based experiment can contain a total of 400 Samples.
- Well Samples are created in a table as part of a 96- or 384-well plate and Tube Samples are created by row.
- Empty Sample spaces are displayed as white squares and Samples that are marked as part of an Experiment show the Experiment and Group color as described in “Experiment/Group labels” on page 216.

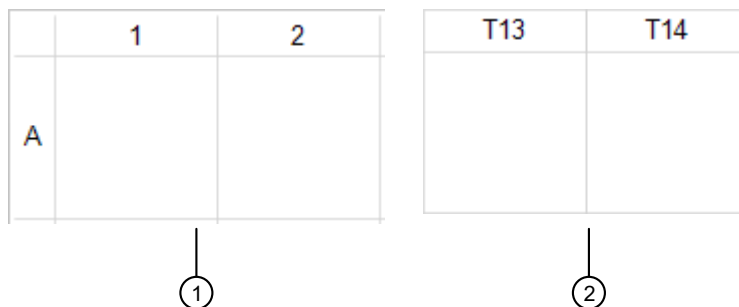


Figure 37 Blank Well and Tube Samples

① Empty Well Sample spaces (A1 and A2)

② Empty Tube Sample spaces (T13 and T14)

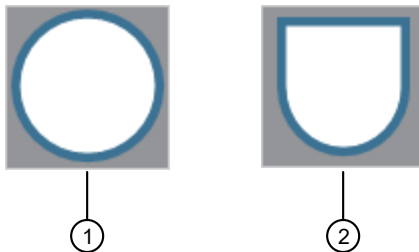


Figure 38 Well and Tube Samples

① Well Sample

② Tube Sample

## Tube Experiment

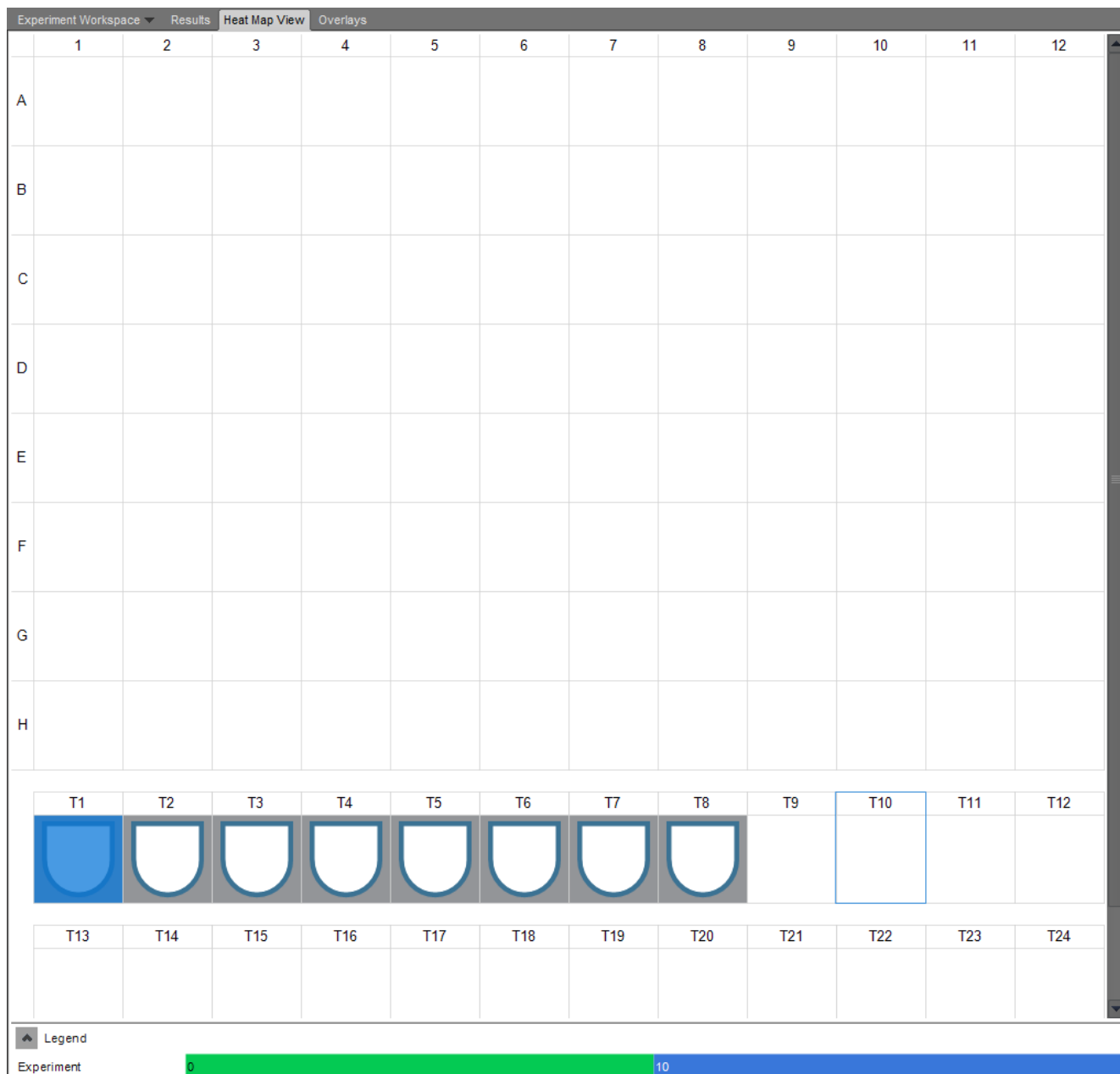
- **Tube Samples** for a new **Tube Experiment** are displayed in rows of 12 until 205 Samples are reached, at which point the view switches to 24 Samples per row.
- Each **Tube Sample** is labeled with its location in the **Heat Map View**, beginning with **T1** and continuing to **T400**, which is the limit to the number of samples a tube- or plate-based experiment can contain.

T1	T2	T3	T4	T5	T6	T7	T8	T9	T10	T11	T12
T13	T14	T15	T16	T17	T18	T19	T20	T21	T22	T23	T24

- The **Heat Map** for a **Tube Experiment** contains the number of **Tube Samples** that were originally created using the **New Experiment dialog** (“New Experiment dialog” on page 606). Any remaining tube spaces in a **Tube Sample** row and the entire row below it are left as empty **Tube Samples**.
- As new **Tube Samples** are generated after the creation of the **Experiment**, the **Heat Map** continues to retain a complete row of empty **Tube Samples** until the total sample limit is reached.

## Plate Experiment

- The **Heat Map** for a new **Plate Experiment** is displayed with an empty table of 96 or 384 wells (depending on the plate type selected), plus the number of **Tube Samples** created in the **New Experiment dialog** (“New Experiment dialog” on page 606).



- 96-well **Plate Experiments** can contain up to 304 **Tube Samples** (for a total of 400 **Samples** per **Experiment**). Any remaining **Samples** are shown as blank **Tube Samples**.
- 384-well **Plate Experiments** can contain up to 16 **Tube Samples** (for a total of 400 **Samples** per **Experiment**). Any remaining **Samples** are shown as blank **Tube Samples**.

- Any **Tube Samples** that are displayed in the **Heat Map** for a **Plate Experiment** are followed by the necessary number of empty **Tube Samples** to complete the row, plus an entire row of empty **Samples**.
- If the **Plate Experiment** contains no **Tube Samples**, then the **Heat Map** shows one complete row of empty **Tube Samples** after the **Well Samples**.
- As new **Tube Samples** are generated after the creation of the **Experiment**, the **Heat Map** continues to retain a complete row of blank **Tube Samples** until the total sample limit is reached.

---

**Note:** **Compensation samples** in tubes are not shown in the **Heat Map**. Therefore, they do not count towards the 400 sample limit.

---

## Navigation and selection

- Double-clicking a **Sample** in the **Heat Map View** loads the **Sample** using the **Experiment-level Workspace**.
- Multiple **Samples** can be selected at one time. The following methods can be used alone or in combination to select the desired **Tube** and/or **Well** samples.
  - Use left-click and drag to lasso a selection of **Tube** and/or **Well Samples**.
  - Click as you hold the **Ctrl** key to select non-contiguous **Tube** and/or **Well Samples**.
  - Click as you hold the **Shift** key to select all wells between the selected points, by row order.
  - Click the **row header** to select that entire row.  
Alternatively, you can use the **Shift+Space** combination to select the entire row, if the row contains at least one selected **Sample**.
  - Click the **column header** to select that entire column.  
Alternatively, you can use the **Ctrl+Space** combination to select the entire column, if the column contains at least one selected **Sample**.
  - Click the top left corner of a **plate diagram** to select the entire plate.  
Alternatively, you can use the **Ctrl+A** combination to select the entire plate, if the plate contains at least one selected **Sample**.

## Set up an Experiment

### Experiment/Group labels

Use the **New Group** (🔍), **New Sample** (🔍), and **Edit Sample** (🔍) buttons to create **Samples** and organize them into **Groups**.

- **Tube** and **Well Samples** have different shape to help identify them more easily.
- Each **Experiment** has a unique identifying color that fills the area surrounding the **Sample** that belongs to it. This color differentiates the **Sample** belonging to that **Experiment** from **Samples** in other **Experiments**.

You can change the **Experiment color** using the **Experiment Information** options in the **Customize panel** (“Experiment information” on page 470) when in the **Heat Map View**.

- Each **Group** has a unique identifying color that marks the inside of the **Experiment color identifier**.

You can change the **Group color** from the **Group Information** options in the **Customize panel** (“Group information” on page 471) when in the **Heat Map View**.

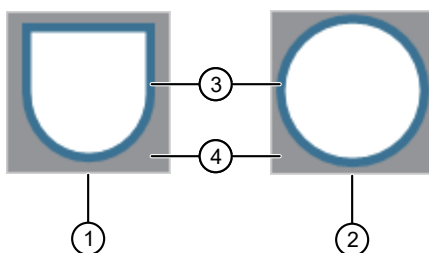


Figure 39 Experiment and Group labels in Heat Map View.

- ① Tube Sample
- ② Well Sample
- ③ Group color
- ④ Experiment color

### Compensation wells

- You can map **Well Samples** as **Compensation Samples** using the **Compensation Setup dialog** as described in “Compensation setup dialog” on page 581.
- **Well Samples** that are mapped as **Compensation Samples** are tagged with a **Compensation Sample icon** **CS** and the name of the parameter they represent. These samples show the **Experiment color** as the **Group color**.

	1	2	3	4	5	6	7
A	CS BL1-A	CS BL2-A	CS BL3-A	RL1-A	RL2-A	RL3-A	UC

## Manual wells

- You can map **Well samples** as **Manual wells** using the **New Manual Well** button in the **Experiment tab** (“Experiment tab controls” on page 95).



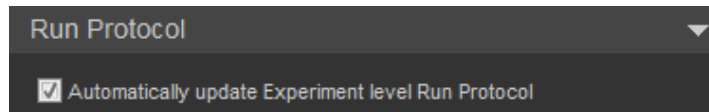
- Well samples** that are mapped as **Manual wells** are tagged with a **manual well icon** **MW**.



## Run Protocol

- You can set up a **Run Protocol** using the **Collection Panel** (Chapter 12, “Collection panel”). Each **Sample** can have a unique **Run Protocol**.

---

**Note:** By default, the **Automatically update Experiment-level Run Protocol** option is selected in the **Run Protocol** in **Collection Panel** (“Automatically update experiment level run protocol” on page 371). You must deselect this option for each sample to have a unique run protocol.



- 
- You can copy and paste a **Run Protocol** using the **Copy Run Protocol**  and **Paste Run Protocol**  buttons in the **Experiment ribbon tab**.  
Alternatively, you can use the hot keys to copy and paste the selected **Run Protocols**.
  - If multiple wells with different **Run Protocol settings** are selected, you can edit individual fields that are in common without altering other settings.
  - Fields with all values in common show the selected value. Fields with differing values are displayed as blank, but are active.

## Legend

The **Legend** for the **Experiment** shows the **Display Mode**, the **color key**, and the **range** that is specified during Heat Map setup (Chapter 14, “Heat map setup panel”).

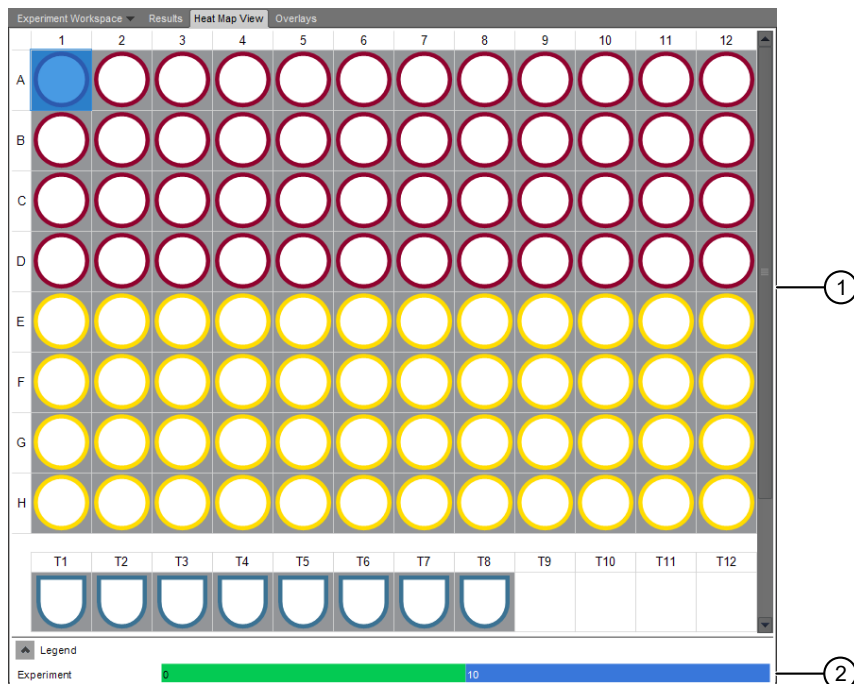


Figure 40 Location of Heat Map View Legend

① Heat Map panel

② Legend (in Threshold mode)

- The **Legend** is attached to the bottom of the **Heat Map panel** and can be minimized, but not closed. It is resized horizontally if the **Heat Map panel** size is changed.
- The **Legend** shows the **Experiment name** in a column on the left, and the **Heat Map** or **Threshold colors** on the right.
- In the **Threshold** mode, **transition values** are shown to the right of the **transition point** (see “Display of results” on page 220).
- In **Heat Map** (i.e., Gradient) mode, **minimum** and **maximum values** are always shown on the **Legend** (see “Display of results” on page 220).
- The value and location of any **transition points** that have been added are displayed to the right of the **transition point** on the **Legend**.



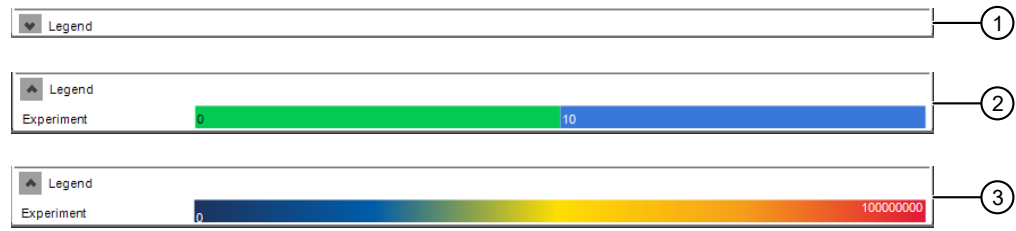


Figure 41 Legend shown in different Display Modes

- ① Legend minimized
- ② Threshold Display Mode
- ③ Heat Map (i.e., Gradient) Display Mode

## Analyze results

### Display of results

Based on the **Heat Map settings** specified in the **Heat Map Setup panel** (Chapter 14, “Heat map setup panel”), each **Sample** with saved data displays a color and, if it is set, the numerical value of the statistic selected for that **Sample**.

#### Threshold display mode

If the **Threshold** is selected for **Display Mode** on the **Heat Map Setup panel**, each **Sample** in the **Heat Map View** displays different color depending on whether the signal from the **Sample** falls above or below the set **threshold value**.

#### Heat Map display mode (i.e., Gradient mode)

If the **Heat Map** (i.e., Gradient mode) is selected as the **Display Mode** on the **Heat Map Setup panel**, each **Sample** in the **Heat Map View** displays the color that corresponds to the appropriate **signal range** set during **Heat Map setup**.

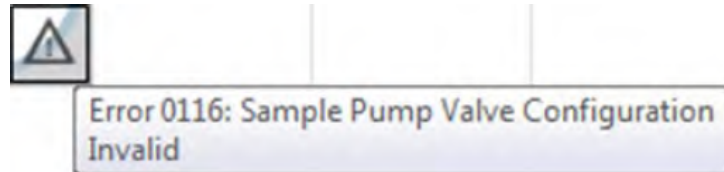
- If the values fall below the **minimum range** set for the sample, the **Heat Map** shows the color representing the **minimum value** and the < symbol.
- If the values exceed the **maximum range** set for the sample, the **Heat Map** shows the color representing the **maximum value** and a > symbol.
- If statistics are selected for display on the **Heat Map**, the < and > symbols are hidden by the displayed statistic.

### Tooltips

- Hover the mouse pointer over a **Well** or a **Tube** to show a **tooltip** with the following information:
  - Gate
  - Statistic
  - Sample
  - Group
  - Time recorded
  - Location
- Hover the mouse pointer over **empty sample spaces** to show a **tooltip** that shows the **Heat Map location** (i.e., plate row and column or tube number).
- Hover the mouse pointer over **Compensation samples** to display a **tooltip** that shows the **compensation parameter** and the **Heat Map location** (i.e., plate row and column or tube number).
- Any fields on the tooltip that are not applicable are left blank.

## Indicators

- If an error causes the Sample run to end prematurely, the **error** icon is displayed in the Heat Map location of the Sample.  
On mouse-over, a tooltip provides more information about the error that was encountered.



- If a bubble is detected during acquisition, the **Bubble detected** icon is displayed in the Heat Map location of the Sample.



- If an error occurs during acquisition, the **generic error** icon is displayed in the Heat Map location of the Sample.





# Sample List view

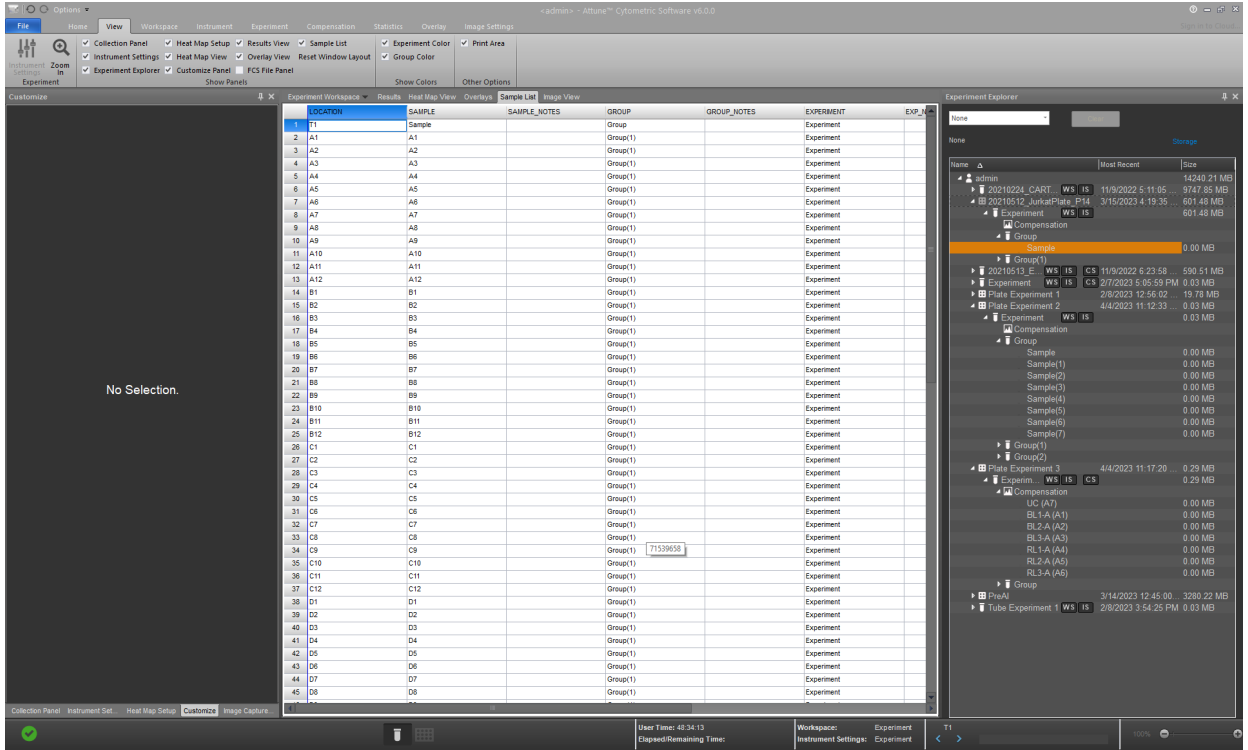
## Overview

The **Sample List** is a grid view of the **Samples** in the current **Experiment**. It displays a table of all **Samples** in the active **Experiment** with columns for the following values:

- Location
- Sample name
- Sample notes
- Group name
- Group notes
- Experiment name
- Experiment notes
- Plate name
- Plate notes
- Plate ID

More columns for **Experiment keyword** are also displayed for all keywords listed in the **Experiment Keywords dialog** (“Experiment Keywords dialog” on page 770).

By default, the **Sample List** is docked to the **Main Application Workspace** (“Main application workspace” on page 56) and is closed. You can open the **Sample List** by selecting it in the **Show Panels** group of the **View** ribbon tab (“Show Panels group” on page 78).



	LOCATION	SAMPLE	SAMPLE_NOTES	GROUP	GROUP_NOTES	EXPERIMENT	EXP_N
1	T1	Sample		Group		Experiment	
2	A1	A1		Group(1)		Experiment	
3	A2	A2		Group(1)		Experiment	
4	A3	A3		Group(1)		Experiment	
5	A4	A4		Group(1)		Experiment	
6	A5	A5		Group(1)		Experiment	
7	A6	A6		Group(1)		Experiment	
8	A7	A7		Group(1)		Experiment	
9	A8	A8		Group(1)		Experiment	

## Sample List properties

### Edit values in the Sample List

- You can edit the following sample information in the **Sample List**:  
Sample name, Sample notes, Group name, Group notes, Experiment name, and Experiment notes
- You cannot edit the Location information (T1, T2, T3... etc; A1, A2, A3... etc).
- Modifications to Group name apply to all Samples in the same Group.
- Modifications to Experiment name apply to all Samples in the Experiment.
- When the **Sample List** for a specific Experiment is in view, extra columns for Experiment keyword are displayed for all keywords listed in the **Experiment Keywords dialog** (“Experiment Keywords dialog” on page 770).  
The title of each column is the name of the keyword.  
When creating a new Experiment, if a default value is specified, that value is populated for the given keyword.  
You can edit the keyword fields for each Sample, or you can leave them blank.
- All changes to the Names and or Notes made in the **Sample List** are propagated throughout the application (i.e., **Customize panel**, **Collection panel**, **Status bar**, **Experiment Explorer**, and the **Statistics** are all updated).
- Name validation occurs when you click elsewhere or when you press the **Enter** key. For the characters and names allowed when editing a field in the **Sample List**, see “Rules for editing the Sample List”.

### Rules for editing the Sample List

- **Name** fields in each grouping must be unique. If names are not unique, a warning message is displayed and the cell reverts to the last unique name used.
- **Sample**, **Group**, **Experiment**, and **Plate names** must be less than or equal to 50 characters long. These fields cannot be blank.
- The **Sample notes**, **Group notes**, and **Experiment notes** can have up to 500 characters. **Note** fields can be left blank.
- Names cannot contain the following illegal characters: / : \* ? " < > | &.
- Names cannot end with a “period” character and cannot be any of the following words: CON, PRN, AUX, CLOCK\$, NUL, COM1, COM2, COM3, COM4, COM5, COM6, COM7, COM8, COM9, LPT1, LPT2, LPT3, LPT4, LPT5, LPT6, LPT7, LPT8, LPT9.
- If a **Sample**, **Group**, **Experiment**, or **Plate name** contains one of these invalid characters or names, the illegal character or name is not accepted and a warning message is displayed. If there is an invalid name, the cell reverts to the last good name used.
- When a column is set to a number type, the cell only accepts digits, a negative character, thousands separator, and the decimal character as per system locale settings. The number of decimal digits allowed is determined by the limit specified in the keyword definition. A number field can be left blank.

---

**Note:** The number fields can only have up to 16 decimals of precision as per the IEEE Standard for Floating-Point Arithmetic (IEEE 754).

---

- If a non-numeric value is entered into a number field, the number field reverts to the last permitted value on validation (i.e., pressing **Enter**, tabbing off the cell, or clicking elsewhere).
- Numeric value formats displayed in the **Sample List** are determined by the system locale settings.

## Navigation in the Sample List with the keyboard

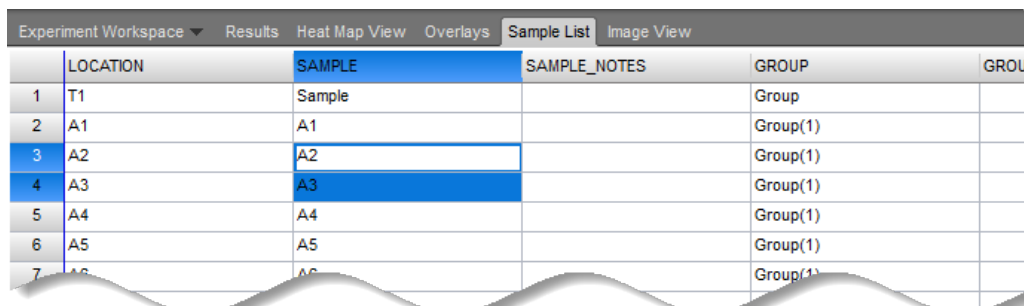
You can navigate the **Sample List** using the keyboard.

- Press the **Tab** key to move from cell to cell from right to left, and to jump to the next row when you reach the last column.
- Press **Enter** to validate the data in the current cell and move to the cell in the next row. When you reach the last row, press **Enter** to jump to the first cell in the next column.
- The **up arrow** and **down arrow** keys move the cell up and down. If the cell is in **Edit mode**, the cell data are validated before advancing to the next well.
- The **left arrow** and **right arrow** keys move the active cell left and right. If the cell is in **Edit mode**, the **left arrow** and **right arrow** keys move the cursor position in the selected text.
- **Page Up** and **Page Down** keys scroll the active view up and down one page at a time.
- To enter a cell into the **Edit mode**, enter any unicode character in that cell. Alternatively, press **F2** to enter the cell into the **Edit mode**.
- You can only edit cells that are in editable columns.

## Selection in the Sample List

Selection in the **Sample List** works similar to selection in Microsoft™ Excel™ application:

- To select a row of cells, click the **row header** or use the **Shift+Space** keyboard combination.
- To select a column of cells, click the **column header** or use the **Ctrl+Space** keyboard combination.
- To select all cells in the **Sample List**, click the **top left corner** of the table or use the **Ctrl+A** keyboard combination.
- Selected cells are highlighted in blue.
- The active cell is shown with a blue border.
- The row and column headers corresponding to selected cells are highlighted with a blue gradient to indicate which rows and columns have selected cells.



	LOCATION	SAMPLE	SAMPLE_NOTES	GROUP	GROL
1	T1	Sample		Group	
2	A1	A1		Group(1)	
3	A2	A2		Group(1)	
4	A3	A3		Group(1)	
5	A4	A4		Group(1)	
6	A5	A5		Group(1)	
7	A6	A6		Group(1)	

## Multiselection in the Sample List

You can select multiple cells in the **Sample List**:

- Click a cell to select it.
- Ctrl+click more cells to select those cells. Focus is set to the last selected cell.
- Shift+click to select the range of cells that are within the rows/columns defined from the first cell to the last cell selected.
- Alternatively, click a cell and drag the mouse to select multiple cells in the **Sample List**. The selected cells are defined by the row/column rectangle created between the start point (last clicked cell) and the cell under the current mouse position. The selection is actively updated as the mouse is dragged.
- To select extra blocks of cells, Ctrl+click, then follow by Shift+click or click and drag.
- You can also select multiple cells using the **keyboard arrows** to select a cell, then using the **Shift** key along with the **left/right/up/down arrows** or **Page Up/Page Down** keys to select the range of cells defined by the row/column rectangle created between the start point (last clicked cell) and the current selected cell.
- If the **Shift** key is not pressed, only a single cell is selected and navigation occurs as described above.
- Press **Ctrl+Shift + an arrow key** to select all cells from the last clicked cell to the limit of the **Sample List** in the direction the arrow key.

## Copy and Paste

You can copy content from the **Sample List** to the clipboard, and then paste it elsewhere in the **Sample List** or to an external spreadsheet such as Microsoft™ Excel™.

Similarly, you can copy content from an Microsoft™ Excel™ spreadsheet, and then paste it into the **Sample List**.

Copy and paste in the **Sample List** works similar to Microsoft™ Excel™ application:

- You can select, copy, and paste multiple columns and rows.
- You can paste only into the editable cells in the **Sample List**. **New Samples** must be added using the **Import Sample List** option in the **Plate** or **Experiment context menus** (“Import Sample List” on page 305 and “Import Sample List” on page 312, respectively).
- To copy the selected cells, use the keyboard shortcut **Ctrl+C** or the **Copy** option in the **Sample List context menu** (“Sample List context menus” on page 227).
- If multiple cells are selected, the destination range must be the same as the source range (i.e., the number of rows and columns must match).
- When pasting contents from clipboard to cells in the **Sample List** that are part of a set (i.e., **Group**, **Group note**, **Experiment**, **Experiment note**, **Plate**, **Plate note**, or **Plate ID**), the items are pasted from top to bottom and left to right.
- The last item pasted in the set is applied to the set.
- If the clipboard contents include invalid characters or illegal names (i.e., system reserved names or names ending in periods), the contents are not pasted into cells where those entries are invalid.



- When pasting into a numeric keyword column and the allowable number of decimal places is exceeded, the number is rounded to the allowable number of decimal places.
- You cannot paste a non-numeric value into a numeric keyword cell.

## Print the Sample List

To print the **Sample List table**, use the print command (**Ctrl+P**) when the **Sample List** is in focus.

Alternatively, select **Print > Sample List** from the **File tab**.

- When printing the **Sample List**, the **Sample List table** is printed from left to right using the current page layout settings (orientation and paper size).
- Columns that do not fit on the same page are printed on the next page.
- Column headers are printed on each page.
- Cell selection does not show up when printed. The column and row header gradients are not printed.

## Sample List context menu

### Sample List context menus

To open the **Sample List context menu**, right-click a cell selected in the **Sample List**.



Figure 42 Sample List context menus – Single and Multi cell

- ① Sample List context menu for a single cell selection
- ② Sample List context menu for multi cell selection

- The **Sample List** context menu has the following options:
- **Edit**: Sets the cell to **Edit mode**. This option is only enabled when the instrument is not acquiring, or the instrument is paused. You can also enter cells into **Edit mode** by double-clicking on them.
- **Clear Contents**: Deletes the contents of the selected cells. This is only enabled for **Sample notes**, **Group notes**, **Experiment notes**, or **Plate notes**, and the instrument is not acquiring, or the instrument is paused.
- **Copy**: Copies the contents of the selected cells to the clipboard. This option is only enabled when the instrument is not acquiring, or the instrument is paused.
- **Paste**: Pastes the clipboard contents to the selected cells. This option is only enabled when there is content in the clipboard to paste, the instrument is not acquiring or paused, and the instrument is not in automation mode.



# Results view

## Overview

The **Results view** displays results from all Samples consolidated into a single table (i.e., **Results table**).

Experiment Workspace															Results															Overlays															Heat Map View														
Drag a column header here to group by that column.																																																											
Group	Workspace	Compensation Source	X Parameter	Gate	Y Parameter	Count	Volume	Concentration	%Total	%Gated	X Mean	Y Mean	X Median	Y Median	X Peak	Y Peak	X SD	Y SD	X %CV	Y %CV	X rSD	Y rSD																																					
Group	Experiment	None	FSC-A	R1	SSC-A	13841	55	251.655	70.448	70.448	438284	226247	436047	214784	4086...	185418	72377.38	77395.62	16.51	34.21	71321.95	7684																																					
Group	Experiment	None	FSC-A	All Events	SSC-A	19647	55	357.218	100...	100.000	458639	264860	438163	219773	4086...	185418	199578...	186772.82	43.52	70.52	112954...	10959																																					
Group	Experiment	None	SSC-A	R2	SSC-H	13795	55	250.818	70.214	99.668	225783	145644	214352	141084	1854...	129914	77050.31	42041.68	34.13	28.87	76201.19	4189																																					
Group	Experiment	None	SSC-A	All Events	SSC-H	19647	55	357.218	100...	100.000	264860	N/A	219773	N/A	1854...	N/A	186772...	N/A	70.52	N/A	109596...																																						
Group	Experiment	None	SSC-A	R1	SSC-H	13841	55	251.655	70.448	70.448	226247	145621	214784	141033	1854...	129913	77395.62	42012.88	34.21	28.85	76643.01	4180																																					
Group	Experiment	None	BL2-A	R3		7142	55	129.855	36.352	51.772	398075	N/A	267276	N/A	3576...	N/A	376090...	N/A	94.48	N/A	355288...																																						
Group	Experiment	None	BL2-A	All Events		19647	55	357.218	100...	100.000	203052	N/A	10094	N/A	-555	N/A	334935...	N/A	164.95	N/A	14213.69																																						
Group	Experiment	None	BL2-A	R2		13795	55	250.818	70.214	70.214	207234	N/A	9586	N/A	-984	N/A	335149...	N/A	161.73	N/A	13027.61																																						
Group	Experiment	None	RL1-A	R4		1	55	0.018	0.005	0.007	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A																																					
Group	Experiment	None	RL1-A	All Events		19647	55	357.218	100...	100.000	1635	N/A	1369	N/A	-33	N/A	1543.37	N/A	94.42	N/A	699.79																																						
Group	Experiment	None	RL1-A	R2		13795	55	250.818	70.214	70.214	1413	N/A	1340	N/A	1333	N/A	534.55	N/A	37.83	N/A	490.74																																						
Group	Experiment	None	RL1-A	-/+	BL2-A	7139	55	129.800	36.336	51.751	1373	398224	1301	267402	1333	357891	509.88	376096.32	37.13	94.44	468.50	35546																																					
Group	Experiment	None	RL1-A	+/-	BL2-A	3	55	0.055	0.015	0.022	4694	44414	4513	22098	4513	10534	323.65	49009.43	6.89	110.35	16.31	1714																																					
Group	Experiment	None	RL1-A	-/-	BL2-A	6644	55	120.800	33.817	48.162	1449	2364	1386	1973	1370	-984	536.61	1748.12	37.03	73.93	506.31	121																																					
Group	Experiment	None	RL1-A	+/-	BL2-A	9	55	0.164	0.046	0.065	4991	3795	4781	3257	4781	1479	819.68	1600.78	16.42	42.18	238.70	133																																					
Group	Experiment	None	RL1-A	All Events	BL2-A	19647	55	357.218	100...	100.000	1635	203052	1369	10094	-33	-555	1543.37	334935.25	94.42	164.95	699.79	1421																																					
Group	Experiment	None	RL1-A	R2	BL2-A	13795	55	250.818	70.214	70.214	1413	207234	1340	9586	1333	-984	534.55	335149.71	37.83	161.73	490.74	1302																																					
Group	Experiment	None	BL1-A	R5		11514	55	209.345	58.604	83.465	445260	N/A	142739	N/A	4333	N/A	470956...	N/A	105.77	N/A	205374...																																						
Group	Experiment	None	BL1-A	All Events		19647	55	357.218	100...	100.000	361489	N/A	39461	N/A	-1464	N/A	455263...	N/A	125.94	N/A	57618.28																																						
Group	Experiment	None	BL1-A	R2		13795	55	250.818	70.214	70.214	371900	N/A	38722	N/A	-2043	N/A	460751...	N/A	123.89	N/A	55311.36																																						
Group	Experiment	None	RL1-A	-/(1)	BL1-A	12175	55	221.364	61.969	88.257	1435	421255	1360	95229	1333	4333	535.44	468823.55	37.31	111.29	496.67	13622																																					
Group	Experiment	None	RL1-A	+/(1)	BL1-A	1	55	0.018	0.005	0.007	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A																																					
Group	Experiment	None	RL1-A	-/(1)	BL1-A	1619	55	29.436	8.240	11.736	1242	970	1201	1761	1246	-2043	476.03	2364.69	38.32	243.89	421.06	91																																					
Group	Experiment	None	RL1-A	+/(1)	BL1-A	0	55	0.000	0.000	0.000	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A																																					
Group	Experiment	None	RL1-A	All Events	BL1-A	19647	55	357.218	100...	100.000	1635	361489	1369	39461	-33	-1464	1543.37	455263.76	94.42	125.94	699.79	5761																																					
Group	Experiment	None	RL1-A	R2	BL1-A	13795	55	250.818	70.214	70.214	1413	371900	1340	38722	1333	-2043	534.55	460751.03	37.83	123.89	490.74	5531																																					
Controls	Experiment	None	FSC-A	R1	SSC-A	10046	99	101.475	64.426	64.426	422597	231820	420499	220605	4045...	205752	81195.27	79532.81	19.21	34.31	88236.94	7979																																					
Controls	Experiment	None	FSC-A	All Events	SSC-A	15593	99	157.505	100...	100.000	394600	239119	397292	208572	4045...	156638	180283...	159317.94	45.69	68.63	156065...	11016																																					
Controls	Experiment	None	SSC-A	R2	SSC-H	9987	99	100.879	64.048	99.413	231039	147811	219970	143270	2057...	130079	78992.54	43165.74	34.19	29.20	79104.12	4430																																					
Controls	Experiment	None	SSC-A	All Events	SSC-H	15593	99	157.505	100...	100.000	239119	N/A	208572	N/A	1566...	N/A	159317...	N/A	66.63	N/A	110164...																																						
Controls	Experiment	None	SSC-A	R1	SSC-H	10046	99	101.475	64.426	64.426	231820	147744	220605	143222	2057...	130079	79532.81	43093.64	34.31	29.17	79797.98	4422																																					
Controls	Experiment	None	BL2-A	R3		3	99	0.030	0.019	0.030	263086	N/A	31736	N/A	18136	N/A	412543...	N/A	156.81	N/A	20163.36																																						
Controls	Experiment	None	BL2-A	All Events		15593	99	157.505	100...	100.000	1165	N/A	879	N/A	-45	N/A	6773.41	N/A	581.33	N/A	502.60																																						
Controls	Experiment	None	BL2-A	R2		9987	99	100.879	64.048	64.048	1044	N/A	869	N/A	1045	N/A	7412.37	N/A	709.80	N/A	398.82																																						
Controls	Experiment	None	RL1-A	R4		193	99	1.949	1.238	1.933	20958	N/A	8144	N/A	6171	N/A	106248...	N/A	506.95	N/A	2235.76																																						
Controls	Experiment	None	RL1-A	All Events		15593	99	157.505	100...	100.000	2994	N/A	1737	N/A	-48	N/A	25802.92	N/A	861.70	N/A	908.83																																						
Controls	Experiment	None	RL1-A	R2		9987	99	100.879	64.048	64.048	2148	N/A	1646	N/A	1654	N/A	14990.60	N/A	697.99	N/A	661.24																																						
Controls	Experiment	None	RL1-A	-/+	BL2-A	2	99	0.020	0.013	0.020	1705	385561	1705	385561	1709	31736	6.36	500383.41	0.37	129.78	6.67	52458																																					
Controls	Experiment	None	RL1-A	+/+	BL2-A	1	99	0.010	0.006	0.010	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A	N/A																																					
Controls	Experiment	None	RL1-A	-/-	BL2-A	9550	99	96.465	61.245	95.624	1696	923	1611	848	1654	964	674.38	415.71	39.76	45.04	619.73	37																																					
Controls	Experiment	None	RL1-A	+/-	BL2-A	434	99	4.384	2.783	4.346	11919	1900	5833	1883	4401	2055	71135.82	633.87	596.82	33.36	1781.34	53																																					
Controls	Experiment	None	RL1-A	All Events	BL2-A	15593	99	157.505	100...	100.000	2994	1165	1737	879	-48	-45	25802.92	6773.41	861.70	581.33	908.83	50																																					
Controls	Experiment	None	RL1-A	R2	BL2-A	9987	99	100.879	64.048	64.048	2148	1044	1646	869	1654	1045	14990.60	7412.37	697.99	709.80	661.24	39																																					
Controls	Experiment	None	BL1-A	R5		15	99	0.152	0.096	0.150	74138	N/A	4077	N/A	4106	N/A	269577...	N/A	363.62	N/A	407.71																																						
Controls	Experiment	None	BL1-A	All Events		15593	99	157.505	100...	100.000	1094	N/A	809	N/A	-25	N/A	8738.21	N/A	798.70	N/A	462.57																																						
Controls	Experiment	None	BL1-A	R2		9987	99	100.879	64.048	64.048	994	N/A	779	N/A	669	N/A	10494.33	N/A	1055...	N/A	345.45																																						
Controls	Experiment	None	RL1-A	-/(1)	BL1-A	20	99	0.202	0.128	0.200	3291	55413	3115	3124	3115	2792	1200.52	233766.29	36.48	421.86	1340.27	43																																					

# Results table

## Results table

The **Results table** shows the statistics associated with gates from the selected Workspaces for **all Samples** in an Experiment.

- **Results table** displays only the gates for the Workspace. Each gate created in the Workspace generates a **results row** in the Results table.
- If **Auto-Refresh** is enabled in the **Home tab**, then results for all files in the Experiment are shown; otherwise, only results for the processed files are shown.

## Primary table

- **Primary table** only displays the statistics selected in the **Statistics tab** (“Statistics tab” on page 103), with each statistic arranged in a separate column in the table.
- By default, columns adjust their width to fit the widest text in the column. The width of a column can be also manually adjusted by dragging the right column divider (i.e., the divider on the right of the column).  
After a column width has been manually adjusted, it does not automatically adjust to fit any new text in the column. Double-clicking the right column divider sets that column back to auto-adjust its width.
- By default, the columns for the primary table are added in the following order:  
Plate, Experiment, Group, Sample, Workspace, Plot Title, Gate, X parameter, Y parameter, Count, Concentration, % Total, % Gated, X mean, Y mean, X median, Y median, X mode, Y mode, X SD, Y SD, X %CV, Y %CV, X rSD, Y rSD, X %rCV, Y %rCV, and Autogate Status.  
Plate, Experiment, Group, and Sample columns display their respective names, and the Workspace column displays the source.
- All columns are added to the table in the order the checkboxes are arranged on the **Statistics tab**, left to right then top to bottom within a Ribbon group. You can reorder the columns by dragging and dropping their headers.
- When a new checkbox is selected, the respective column is inserted in the correct position in the table.
- You can remove the columns from the table by deselecting them in the **Statistics tab** or by dragging a column header outside of the column header area. When dragging a column header, an **X** indicates that it will be removed.
- If a statistical value cannot be calculated, **N/A** is displayed instead of a numerical value.
- If a value is not applicable to the row (e.g., Y values for a histogram gate), it is left blank.
- The table can be sorted by any column in an ascending or descending order (case insensitive) by clicking the column title.

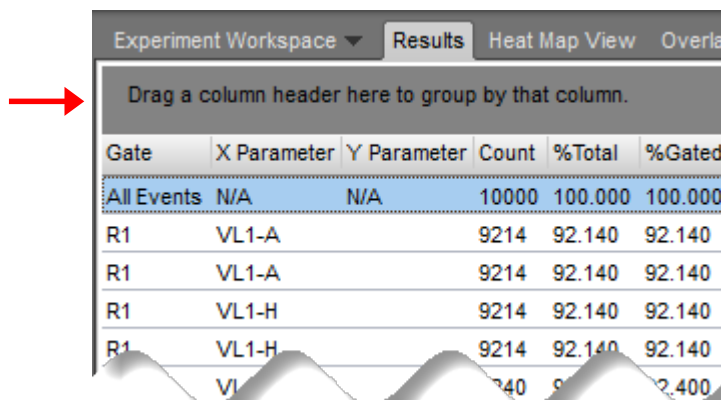
## Select rows

- Left or right mouse click a **Results row** to select the entire row.
- Hold down the **Ctrl** or **Shift** keys to select multiple rows.
- **Ctrl+A** combination selects all the rows in the table.
- **Ctrl+Shift+End** combination selects all rows from the currently selected row to the end of the table.
- **Ctrl+Shift+Home** combination selects all rows from the currently selected row to the top of the table.

## Group results

### Group results by column headers

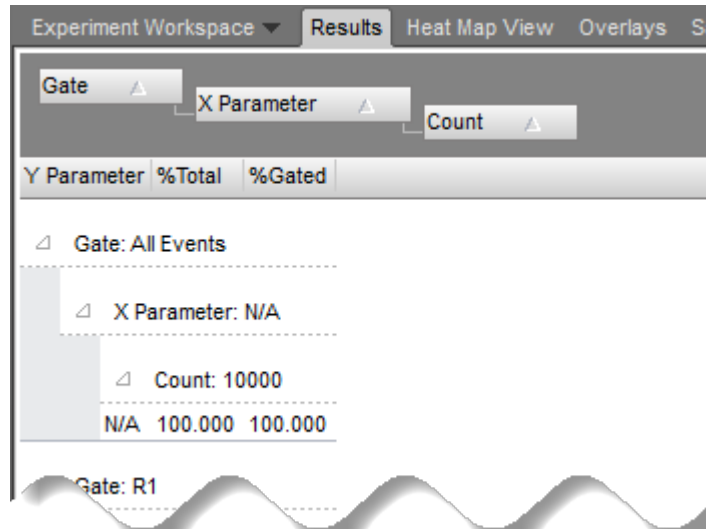
- Results can be grouped by the values displayed in the column headers by dragging a column header from the **Results table** onto the area above the table that is labeled **Drag a column header here to group by that column.**






The screenshot shows the 'Results' view in the software. At the top, there are tabs for 'Experiment Workspace', 'Results', 'Heat Map View', and 'Overlay'. Below the tabs is a grey instruction box that says 'Drag a column header here to group by that column.' A red arrow points to this instruction box. Below the instruction box is a table with the following data:

Gate	X Parameter	Y Parameter	Count	%Total	%Gated
All Events	N/A	N/A	10000	100.000	100.000
R1	VL1-A		9214	92.140	92.140
R1	VL1-A		9214	92.140	92.140
R1	VL1-H		9214	92.140	92.140
R1	VL1-H		9214	92.140	92.140
VI			9214	92.140	92.400

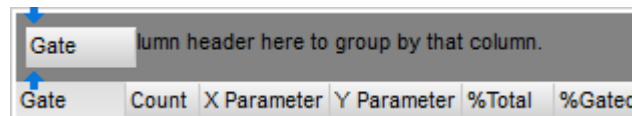
- You can place more than one header in the **Drag a column header here...** area, which groups the results by all the headers placed there. The headers are linked by arrows denoting the sort order:



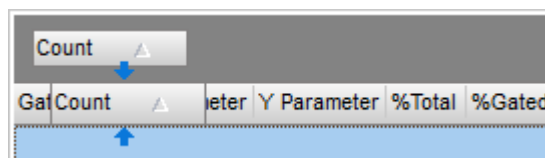
- You can change the sort order by dragging a header in a different position on the group control.
- Clicking on a header arrow (  ) toggles the sorting order between ascending and descending in the column.
- To collapse the information displayed in the groups, click the  button; to expand it, click the  button.

## Add and remove column headers

- When a **header** is dragged, two **blue divider arrows** indicate where the **column header** will be placed in the **Drag a column header...** group area.



- To remove a **header** from the grouping, drag it back to its position in the **Results table**. When a **column header** is dragged back to the **Results table**, two **blue divider arrows** indicate where the header will be placed.



Alternatively, drag the header off the table to remove it from the **Drag a column header...** group area.

## Moving versus copying headers

- Dragging a header into the **Drag a column header...** group area removes the **header** from the **Results table** and places it in the **group area**.  
Dragging the **header** back to the **Results table** removes it from the **group area** and places it back in the **Results table**.
- If you hold down the **Ctrl** key as you drag a **header**, the **header** is copied instead of moved.

## Examples of grouped results

Gate						
Group	Sample	Count	Volume	Concentration	%Total	%Gated
△ Gate: All Events						
Group	Spleen Control	2239344	212.0	10562.9	100.0	100.0
Group	Spleen Control	2239344	212.0	10562.9	100.0	100.0
Group	Spleen Control	2239344	212.0	10562.9	100.0	100.0
Group	Spleen Control	2239344	212.0	10562.9	100.0	100.0
Group	Spleen Control	2239344	212.0	10562.9	100.0	100.0
Group	Spleen Control	2239344	212.0	10562.9	100.0	100.0
Group	Spleen Control	2239344	212.0	10562.9	100.0	100.0
Group	Spleen Control	2239344	212.0	10562.9	100.0	100.0
Group	Spleen Control	2239344	212.0	10562.9	100.0	100.0
△ Gate: CD3+						
Group	Spleen Control	500075	212.0	2358.8	22.3	43.1
Group	Spleen Control	500075	212.0	2358.8	22.3	22.3

Figure 43 Example of results grouped by Gate

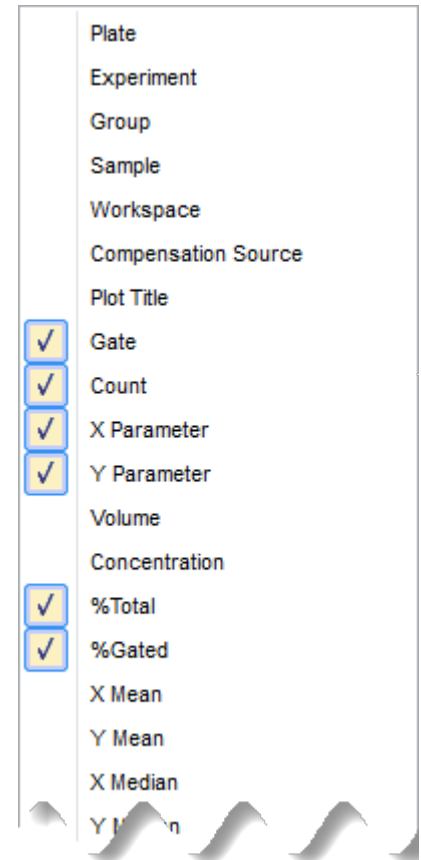
Experiment Workspace ▾ Results Overlays Sample List Heat Map View						
Gate						
X Parameter						
G.	Sample	Count	Volume	Concentration	%Total	%Gated
△ Gate: All Events						
△ X Parameter: BL1 - FJComp-BL1-A						
G..	Spleen Control	2239344	212.0	10562.9	100.0	100.0
△ X Parameter: FSC - FSC-A						
G..	Spleen Control	2239344	212.0	10562.9	100.0	100.0
△ X Parameter: FSC - FSC-H						
G..	Spleen Control	2239344	212.0	10562.9	100.0	100.0

Figure 44 Example of results grouped by Gate and X Parameter

## Results table column header context menu

Right-click a **column header** in the **Results table** to open the **Results table column header context menu**.

- The **column header context menu** contains all the **statistics** that can possibly be displayed in the **Results table**.
- The **statistics** that are already present in the **Results table** are checked in the context menu.
- To add new columns to the **Results table**, check the corresponding item in the **column header context menu**.
- To remove columns from the **Results table**, deselect the corresponding item in the **column header context menu**.
- The **column header context menu** can be displayed when there are no columns in the table or when you click a blank area of the header.



## Overview

**Overlays** provide a graphical method for comparing Experiment data, in which selected plots are superimposed for a direct visual comparison of overlapping data.

- Each **Overlay** consists of an Overlay plot (“Overlay plots” on page 237) and an associated **Overlay gallery** containing the **Source plots** for the **Overlay** (i.e., **Gallery plots**; see “Gallery plots” on page 241).
- **Overlays** are viewed and modified in the **Overlays tab**.

## Overlays tab

**Overlays** tab is displayed in the **Application Area** on the **Main Application Workspace** (“Main application workspace” on page 56) and enables direct comparison of **Histogram** or **dual-parameter plots**.

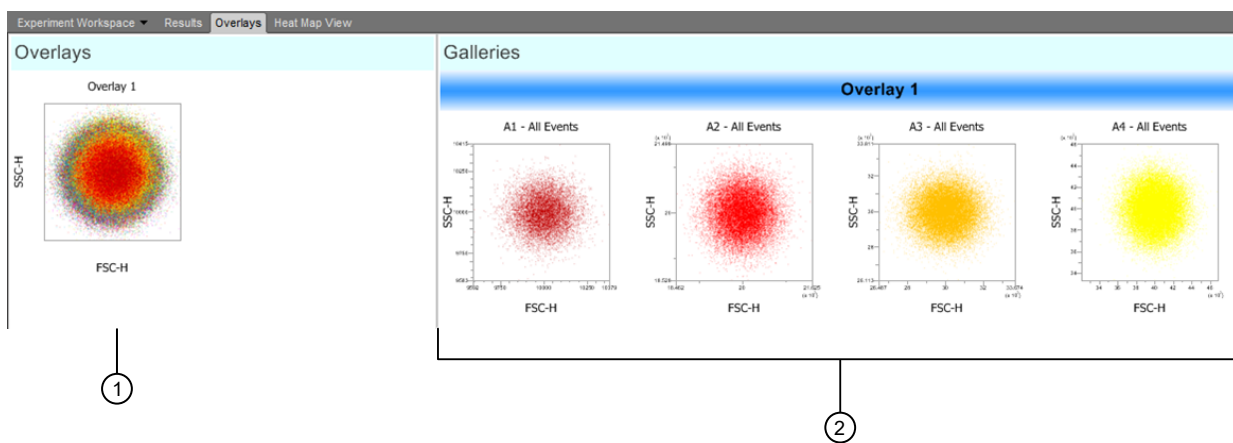


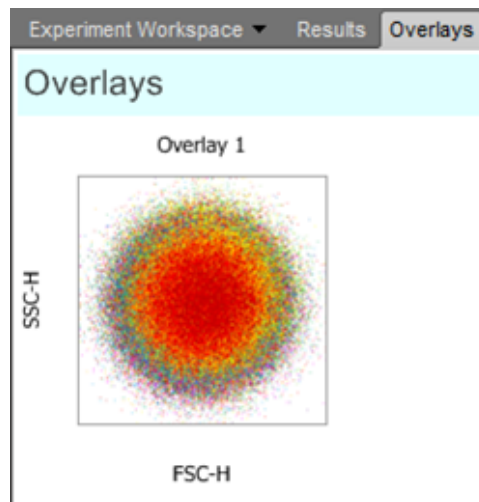
Figure 45 Overlays tab

- ① Overlays area
- ② Galleries area
- The **Overlays** tab is divided vertically into two windows by a splitter bar.
  - The left window of **Overlays** is the **Overlays area** (“Overlays area” on page 235)
  - The right window is the **Galleries area** (“Galleries area” on page 236)
- The docking behavior of the **Overlays** tab is described on “Overlay tab” on page 110.
- There is no limit to the number of **Overlays** that can be added to the **Overlays** tab.
- When the **Customize Overlay panel** (“Customize Overlay options” on page 474) is visible, the panel contents are updated based on the selected plots. The customize panel enables you to customize the appearance of **Overlay** and **Gallery plots**.



## Overlays area

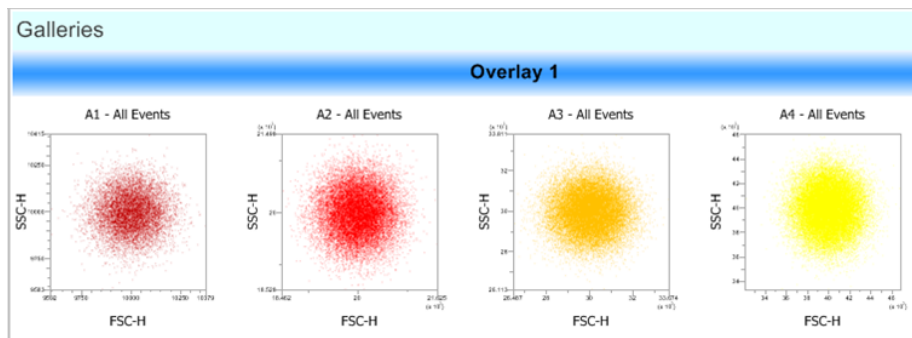
The **Overlays area** displays the **Overlay plots** (“Overlay plots” on page 237).



- To resize the **Overlays area** move the **splitter bar** to the left or to the right.
- **Plots** in the **Overlays area** automatically wrap, when needed. By default, the divider is set so that only one **Overlay plot** is shown per row.  
When needed, **vertical** and **horizontal scroll bars** become available on the **Overlays area** to enable the viewing of all existing plots.
- When new plots are added to an existing **Overlay** or if a new **Overlay** is created by sending plots from the **Workspace**, the most recently created and/or modified **Overlays** are selected when viewing the **Overlays tab**.
- When one or more **Overlay plots** are selected, the titles for the corresponding galleries are highlighted in blue.
- Overlay plots are displayed 240 × 240 pixels. If **legends** are enabled, the **Overlay plots** are resized to 240 × 480 pixels (H × W).

## Galleries area

The **Galleries area** contains the **Galleries** of **Source plots** (i.e., **Gallery plots**; see “Gallery plots” on page 241) for each Overlay plot displayed in the Overlays area.



- Each **Gallery** starts on a new line and has the same title as the **Overlay** whose **Source plots** it contains. The **Gallery title** is shown above the plots and is always visible, even when the plots are horizontally scrolled.
- **Horizontal** and **vertical scroll bars** are available, when needed.
- The **Galleries area** can be resized by moving the **splitter bar** to the left or to the right.
- Selecting any plot in a Gallery highlights the **Gallery title** in blue.
- To resize the Plots in the **Galleries area**, use the **Size Slider** in the lower right corner of the on the **Application Status Bar** (“Size slider” on page 70) or hold down **Ctrl** on the keyboard and turn the **mouse scroll wheel**.
- **Gallery plots** are displayed in 240 × 240 pixels. If **legends** are enabled, the **Gallery plots** are resized to 240 × 480 pixels (H × W).

---

**Note:** Galleries view cannot be printed.

---

## Overlay plots

**Overlay plots** enable the visualization and comparison of overlapping data by superimposing the selected Histogram or of dual-parameter plots. The **Overlay plots** are displayed in the **Overlays area** (“Overlays area” on page 235).

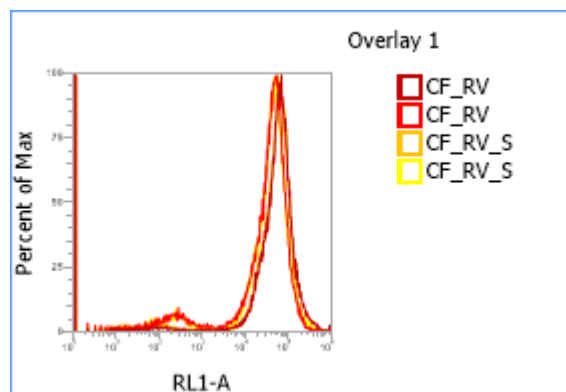
### Overlay plot modes

Overlays plots have two display modes: **Overlay mode** and **3D mode**.

To switch between **Overlay** and **3D modes**, select or deselect the **View 3D** option in the **Customize Overlay panel** (“Customize Overlay options” on page 474). The selected display mode only applies to the currently selected **Overlay**.

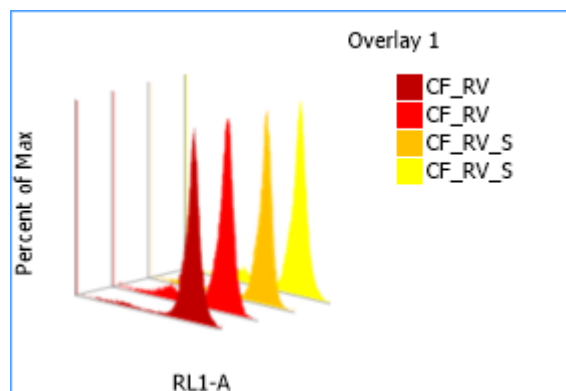
#### Overlay mode

- The **Overlay mode** is the default display mode, where the **Overlay plots** are shown superimposed over one another.

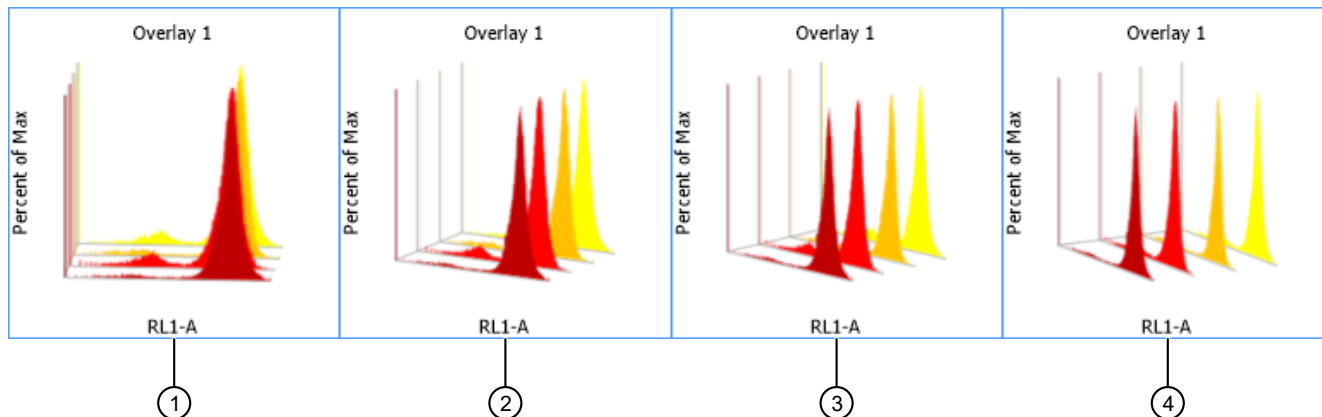


#### 3D mode

- In the **3D mode**, plots are displayed three dimensionally at a selected angle and are separated by a set distance. The distance between the plots is calculated based on the number of plots being overlaid.



- In this mode, you can rotate the plot by selecting the desired angle from the **Angle (deg)** dropdown control in the **Customize Overlay panel** (“Angle (deg)” on page 480).



- ① Angle at 5 degrees
- ② Angle at 30 degrees

- ③ Angle at 45 degrees
- ④ Angle at 60 degrees

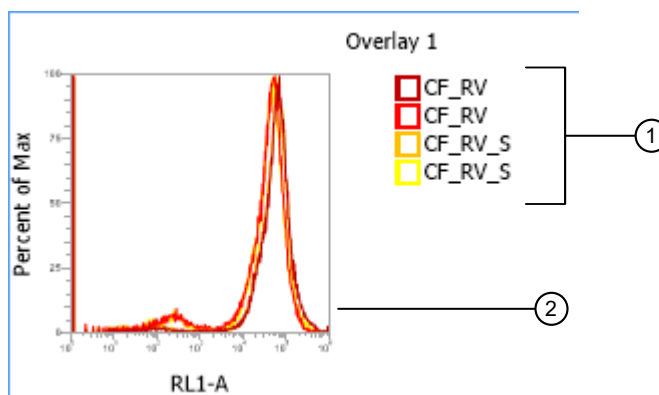
## Overlay plot characteristics

### Plot title

- Overlay plot titles** are defined using the **Text** options on the **Customize Overlay panel** (“Customize Overlay options” on page 474).

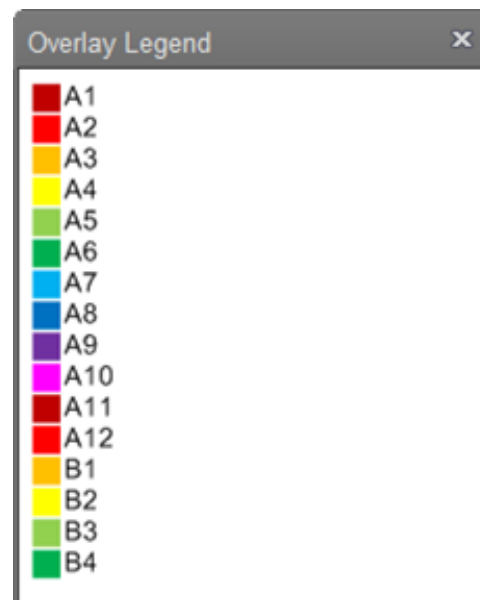
### Legend

- By default, the **legend** is not displayed. However, you can select to display the **legend** by selecting the **Show Legend** option in the **Customize Overlay panel** (see “Show legend” on page 475). The option to display the legend can also be set using the **Set overlay legend** option in the **Overlay Builder** as described in “Show overlay legend” on page 268).
- By default, the **legend** shows the name of each **Sample** in the **Overlay**. The **Parameter name** can also select to show the legend with or without the **Sample name** using the **Customize Overlay panel** (“Show legend” on page 475).



- ① Legend
- ② Overlay plot

- The color associated with each sample in the legend corresponds to the data color of that sample on each plot.  
Name labels of dual parameter plots in the legend are represented by solid colored rectangles. Name labels of single parameter plots in the legend are represented by unfilled rectangles with colored borders unless the single parameter plot is itself filled. In that case, the legend is also displayed as filled.
- If all the plots in the **Overlay** cannot be represented on the **legend** due to insufficient space, the **See All hyperlink** is shown below the last legend item shown.
- Click the **See All hyperlink** to open a floating **Overlay Legend list** that shows all the plots on the **Overlay**.
- Click the **X** (close) button on the floating list to close it.



## Axis labels

- In the **Overlay mode**, an **axis label** is shown only if it is the same for all the plots in the **Overlay**. If an **Overlay** is created from plots where a parameter is different for one of the axes, no **axis label** is shown for that axis.
- To edit the **axis labels** on the **Overlay plot**, use the **Text** options in the **Customize Overlay panel** ("Text" on page 483).

## Axis tick marks

- For the selected **Overlay plots**, you can specify the presence of **X-axis** and/or **Y-axis tick marks** by selecting the **Show X-Axis Ticks** and **Show Y-Axis Ticks** options in the **Customize Overlay panel** (see "Show X-Axis and show Y-Axis ticks" on page 476).
- By default, both the X-axis and Y-axis tick mark display options are enabled.
- Changing tick mark display options for an **Overlay plot** updates all the plots in the associated gallery.
- By default, the tick marks are displayed on **Overlay plots** when all the **Source plots** have the same scale and range. If the scale ranges are different, the tick marks are not displayed, even if they are enabled on the **Customize Overlay panel**.

## Gates

- To add **Gates** to **Overlay plots** in **Overlay mode**, use the **Gating Tools** on the **Overlay ribbon tab** (“Overlay tab” on page 110).
- When a gate is drawn on an **Overlay plot**, the same gate is drawn on all the plots in the associated **Overlay gallery**.

## Re-ordering plots

- To reorder **Overlay plots** in the **Overlay mode**, drag-and-drop the selected plots.  
When you reorder **Overlay plots**, their respective **Overlay galleries** are also reordered to match the order of the **Overlay plots**.
- When dragging plots, a blue vertical line indicates the insertion location of the plots.
- When you select and drag multiple plots, they are grouped together and can be moved and inserted as a group.
- If more than 3 plots are selected, a plus (+) icon is shown next to the plot images indicating the presence of added plots being dragged.
- You can move plots only when and where the **insert indicator** appears; otherwise the plots return to their original positions when released.

## Gallery plots

**Gallery plots** are the **Source plots** that are overlaid to generate the **Overlay plot**. **Gallery plots** are shown in the **Galleries area** (“Galleries area” on page 236).

### Gallery plot characteristics

#### Plot title

- **Gallery plot titles** are defined using the **Text** options on the **Customize Overlay panel** (“Text” on page 483).

#### Axis labels

- **Gallery plot axis labels** can be edited using the **Text** options on the **Customize Overlay panel** (“Text” on page 483).

#### Axis tick marks

- By default, both the X-axis and Y-axis tick mark display options are enabled.
- Tick mark visibility for the selected **Gallery plots** can be changed using the **Show X Axis Ticks** and **Show Y Axis Ticks** checkboxes on the **Customize Overlay panel** (“Show X-Axis and show Y-Axis ticks” on page 476).

#### Order of plots in the Galleries area

- The order of the plots in the **Galleries area** (left to right) determines the z-order (front to back) of plots in the corresponding **Overlay plot**.
- Plots can be rearranged in each **Overlay gallery** by dragging and dropping. When plots are rearranged in an **Overlay gallery**, they are rearranged in the respective **Overlay plot** as well.
- When dragging plots, a vertical blue line highlights the insertion location of the plots.
- When multiple plots are selected and dragged, they are grouped together, and moved and inserted as a group.
- If more than 3 plots are selected, a plus (+) icon is shown next to the plot images indicating the presence of added plots being dragged.
- Plots can only be moved when and where the **insert indicator** appears; otherwise they return to their original positions when released.

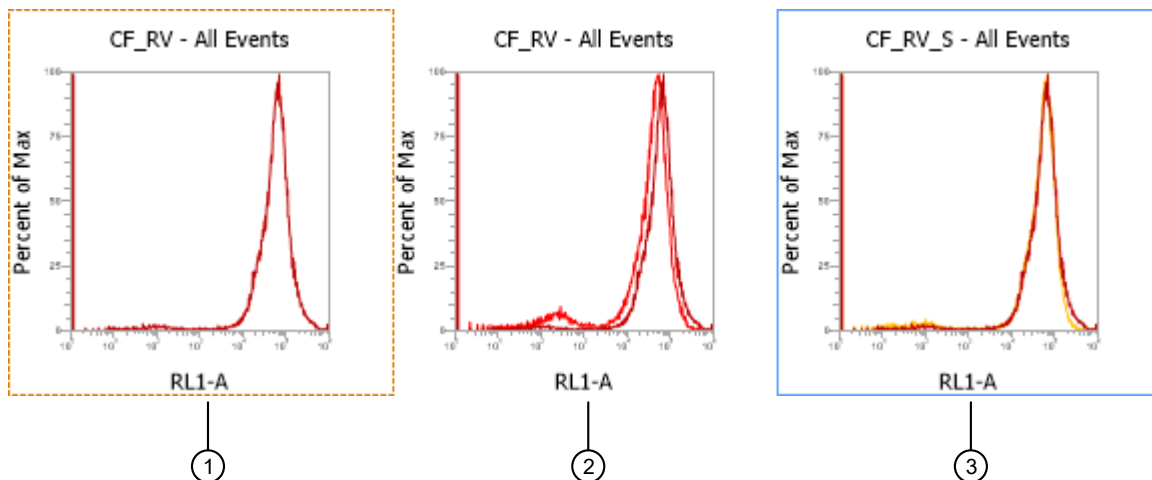
---

**Note:** **Gallery plots** cannot be printed.

---

### Control plot

- In the **Galleries area**, one **Gallery plot** for each single parameter overlay can be selected to be a **Control plot**.
- The **Control plot** is shown as an overlay on all single parameter plots present for a specific overlay in the gallery. The specified **Control plot** is indicated with an orange dashed line around it.



- ① Control plot
- ② Gallery plot
- ③ Gallery plot (selected)

- By default, a **Control plot** is not defined.
- The **Control plot** is selected by clicking the **Set as Control** button in the **Customize Overlay panel** (“Set as Control” on page 481).
- If there is already a **Control plot**, defining a different plot as a new **Control plot** replaces the previous plot as the control.
- When the **Move to Front** option in the **Customize Overlay panel** (“Set as Control” on page 481) is selected, the **Control plot** is shown overlaid in front of the **Test plot** on each **Gallery plot**.  
By default, the **Move to Front** option is not selected and the **Test plots** are shown in front of the **Control plot**.
- When a **Control plot** has been defined, the software enables comparison statistics to be calculated (see “Overlay calculations” on page 243).



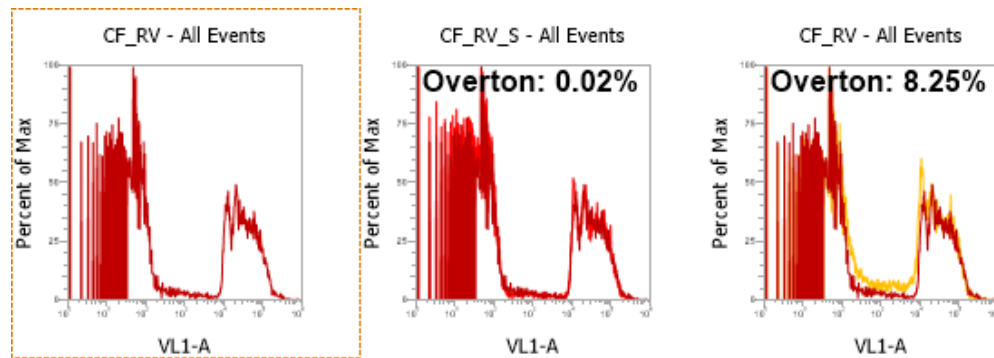
## Overlay calculations

- When a **Control plot** has been defined, the data sets can be compared statistically through selection of the **Calculation method** options in the **Customize Overlay panel** (“Calculation method options” on page 482).

The available options are **None**, **Overton** (“Overton’s cumulative statistics” on page 937) and **KS** (Kolmogorov-Smirnov) (“Kolmogorov-Smirnov (K-S) method” on page 243). Both **Overton** and **KS** (Kolmogorov-Smirnov) methods are used to determine the statistical difference between two samples.

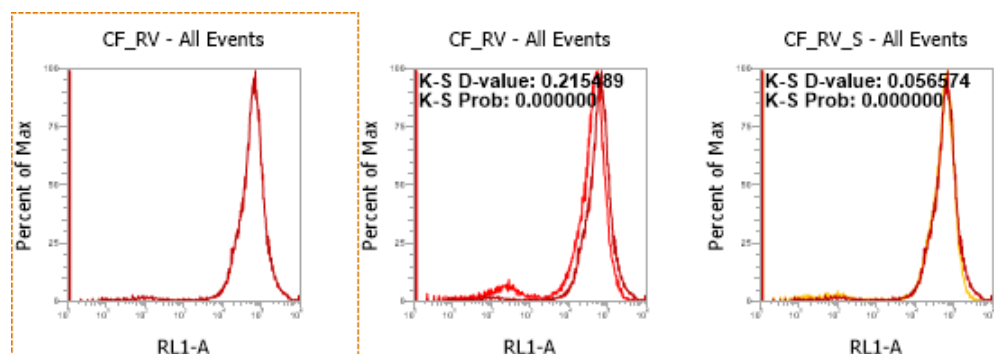
### Overton method

- Overton** option compares the **Control plot** and each **Gallery plot** using the **Overton’s cumulative statistics calculation** (“Overton’s cumulative statistics” on page 937).
- The resulting value indicates the percent difference between the **Test sample** and the **Control sample** and is shown on each test plot. The calculations are not shown on the **Control plot**.



### Kolmogorov-Smirnov (K-S) method

- KS** option compares the **Control plot** and **Test plot** using the Kolmogorov-Smirnov test (“Kolmogorov-Smirnov test” on page 937) to determine the probability that the two plots are the same. Results are displayed on the plot.



- The lower the calculated **D-value**, the more likely that the plots are the same. The lowest possible **D-value** is 0, in which case the plots are identical.
- The **Probability value** is a value between 0 and 1. The greater the **Probability value**, the greater is the probability that the plots are the same.

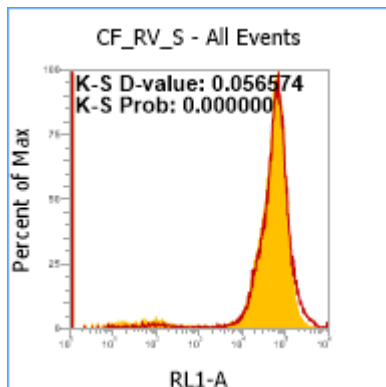
---

**Note:** For more information about the **Overton cumulative statistics calculation** and the **Kolmogorov-Smirnov (K-S) test**, see Appendix C, “Data management in Attune™ Cytometric Software”).

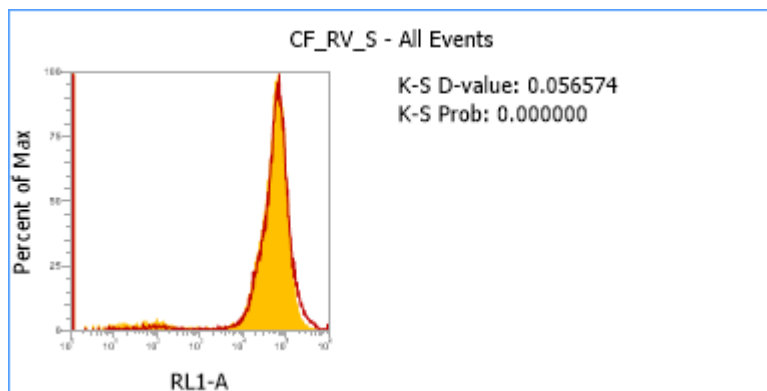
---

## Display of statistics

- **Overlay calculation statistics** (“Overlay calculations” on page 243) and **Overlay gate statistics** (“Customize Overlay options” on page 474) can be shown on the plot or to the side of the plot. By default, they are shown on the plot.



- When the **Show Legend** option (“Show legend” on page 475) is selected on the **Customize panel**, statistics are displayed on the side of the plot as a legend.



- If all the plots and statistics cannot be shown on the legend due to insufficient space, a hyperlink titled **See All** is displayed below the last legend item shown. Clicking the **See All** hyperlink displays a floating list that shows all the plots on the **Overlay**. Clicking the close button (**X**) on the floating list closes it.
- Overton statistics always show two decimal places on the plot. All other statistics show decimal places as specified in the **Options dialog** (“Stats Options” on page 665).

## Gates

- **Gates** are created on a selected plot using the **Gating Tools** available on the **Workspace ribbon tab** (“Gating Tools group” on page 83).
- **Gates** created on a **Gallery plot** are displayed on all the plots in the **Overlay gallery** and on the related **Overlay plot**, if the **Overlay plot** is in **Overlay** mode and not in **3D** mode (“Overlay plots” on page 237).
- **Gates** are not displayed on **Overlay plots** that are in the **3D** mode (“Overlay plots” on page 237).
- **Gates** created on an **Overlay plot** in the **Overlay** mode are reflected on all the plots in its **Gallery**.
- The maximum number of gates for an Overlay is 5.  
If you attempt to add a gate after the maximum number of gates on that plot has already been reached, a **no entry** symbol (⊘) is displayed, indicating that the extra gate cannot be created.
- Only one **Quadrant gate** can be drawn per Overlay. A Quadrant counts as a single gate, therefore allowing 4 extra, non-Quadrant gates on the same plot.
- **Autogates** are not supported in **Overlay plots**.
- **Gates** are drawn in black with no fill and are assigned default names in the order of their creation.
- When a gate is moved or resized on any plot in the **Overlay**, it is updated accordingly in all the remainder of the plots in the same **Overlay**.
- **Statistics** are shown on the **Gallery plot** as described in “Display of statistics” on page 244; they are not shown on the **Overlay plot**.
- The default statistic displayed is **%Gated** (“Event Statistics group” on page 108). You can change the statistics displayed by selecting the desired one from the **Statistics ribbon tab** (“Statistics tab” on page 103).
- The decimal places for the statistics are as set in the **Options dialog** (“Stats Options” on page 665). The statistics shown persist on a per Experiment basis.
- If the statistic cannot be calculated, **N/A** is displayed.
- **Gates** on the **Overlay tab** do not affect the **Source plots** on the **Workspace** and do not count towards the maximum numbers of gates.

# Overlay properties

## Overlay settings

### Overlay title

- The **Overlay title** is unique for each **Overlay**, and it distinguishes the **Overlay plot** and its respective **Overlay gallery** from other existing Overlays.
- The default title for a new Overlay is **Overlay n**, where **n** is a numerical suffix (first available integer) to ensure that the title is unique. If an **Overlay** is deleted or renamed, the number designation is reused.

### Overlay plot colors

- By default, all dual-parameter plots (i.e., **Dot**, **Density**, and **Precedence Density plots**) in an **Overlay** are displayed in a solid color.
- By default, all single-parameter plots are displayed as a solid line, without fill.
- The color of a plot in an **Overlay** is determined by the defaults specified in the **Color and Themes** tab of the **Options dialog** (“Colors and Themes” on page 647).  
Each new plot in the **Overlay** is assigned the next available color listed in the **Gates and Sample Overlays** group in the **Options dialog** ▶ **Color and Themes** tab.

### Range

You cannot change the range of individual plots on the **Overlays tab**.

### Parameter

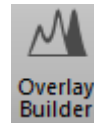
You cannot change parameters of plots on the **Overlays tab**.

### Persistence of Overlay settings

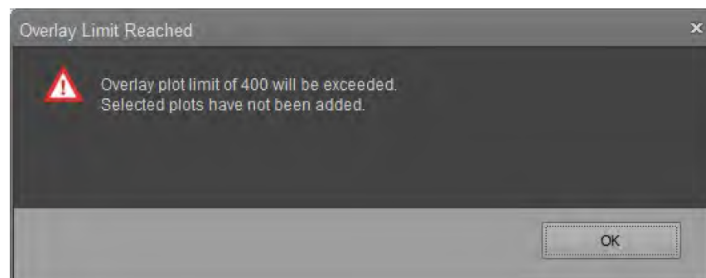
All plots, options, and settings defined on the **Overlays tab** persist at the Experiment level.

## Add plots to Overlays

- To send a **Source plot** from a **Workspace** to a new or existing **Overlay**:
  - Right-click any empty area (i.e., white space) in the **Workspace** to open the **Workspace view context menu**.  
From the **Workspace view context menu**, select **Send All Plots to Overlay** (“Workspace view context menu” on page 185) to overlay all plots in the **Workspace** of the single **Sample** on top of each other.
  - Right-click any empty area within the boundary of a single **Plot** to open the **Plot context menu**.  
From the **Plot context menu**, select **Send to overlay** (“Plot context menu” on page 189), then select the appropriate **Overlay** from the dropdown menu. The selected plot is added to the selected **Overlay**.
  - Alternatively, click the **Overlay Builder** button on the **Home ribbon tab** or the **Overlay ribbon tab** to create an **Overlay** using the **Overlay Builder dialog** (“Overlay Builder” on page 260).



- Each **plot** is added to the selected **Overlay plot** and to its respective **Overlay gallery**.
- By default, **plots** added to a new **Overlay** are arranged in the order of addition.  
When new **plots** are added to an existing **Overlay**, they are added to the end of the **Gallery**.
- By default, new **Overlay plots** are displayed in the **Overlay mode** (“Overlay plots” on page 237).
- **Gates** present on the **Source plots** in the **Workspace** are not added to the **Overlay**.
- **Range** and **Scale** settings of the **Source plots** are applied to the **Overlay plots**. The **Scale** setting of the newly added **Source plot** always takes precedence over any existing scaling set in the **Overlay**.
- An **Overlay** can contain a maximum of 400 **Source plots**. If the addition of new plots causes the **Overlay** to exceed the 400 plot limit, the software displays a warning dialog.



Clicking **OK** closes the warning dialog without adding the new plots to the **Overlay**.

## Compare Plots with different Scale and Range settings

- **Plots** with different **Scale** settings (i.e., **Linear**, **Logarithmic**, or **Hyperlog™**) or different scales of the same type can be compared on the same **Overlay**.

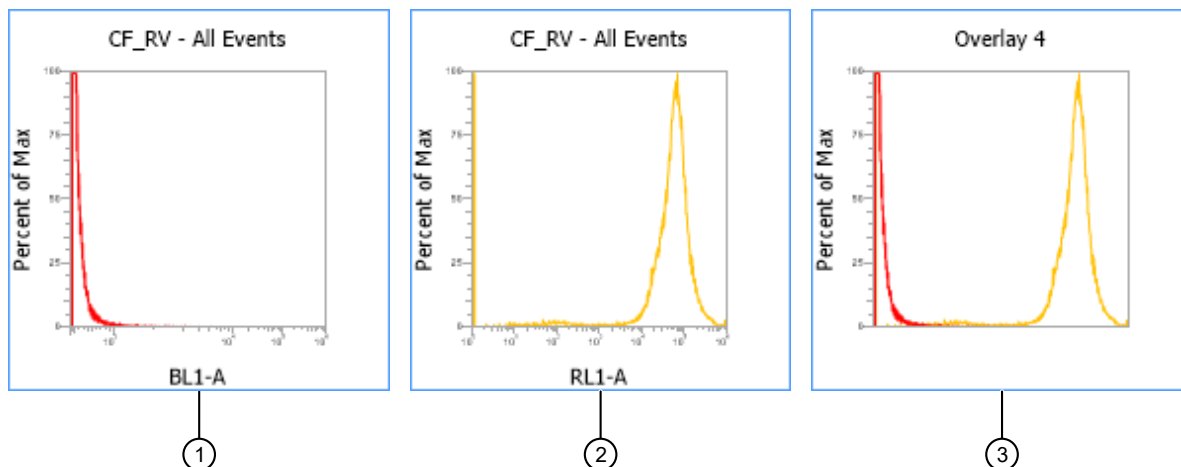


Figure 46 Overlay of plots with different Scale and Range settings

- ① Plot with the x-axis in Logarithmic scale
  - ② Plot with the x-axis in Hyperlog™ scale
  - ③ Overlay of the two plots (Plot 1 and Plot 2) with different x-axis scales
- **Plots** with different **Range** settings can be compared on the same **Overlay**. When comparing plots with different **Range** settings, **axis labels** behave as described in “Add plots to Overlays” on page 247.

---

**Note:** When comparing plots with different scale and range, the settings of individual plots remain unchanged.

---

## Select plots

### Select a single plot

- To select a single **Overlay plot** or **Gallery plot**, click the desired plot. The selected plot is surrounded by a **light blue rectangle**.

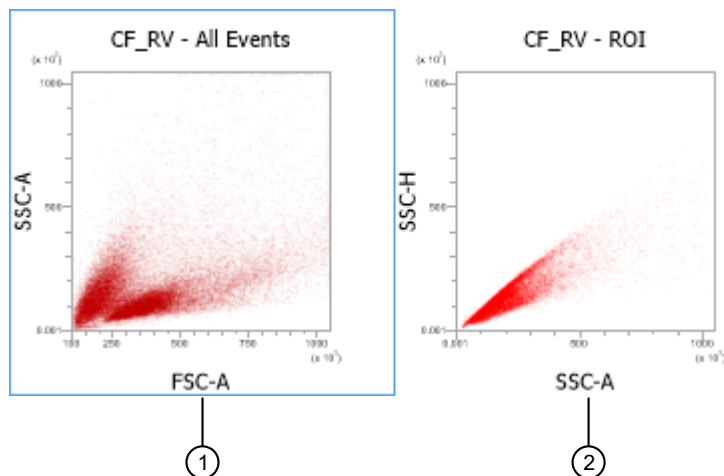


Figure 47 Selection of a single plot

- ① Selected plot surrounded by a light blue rectangle
- ② Plot that is not selected

### Select multiple plots

- To select multiple plots, click the desired plots as you hold down the **Ctrl** key. Objects which are part of the multiselect are bounded by **dark blue rectangles**.

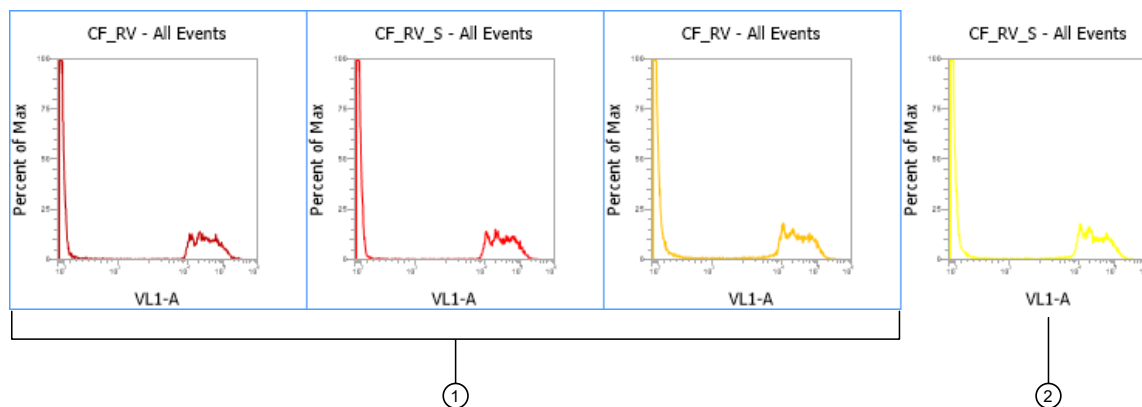


Figure 48 Selection of multiple plots

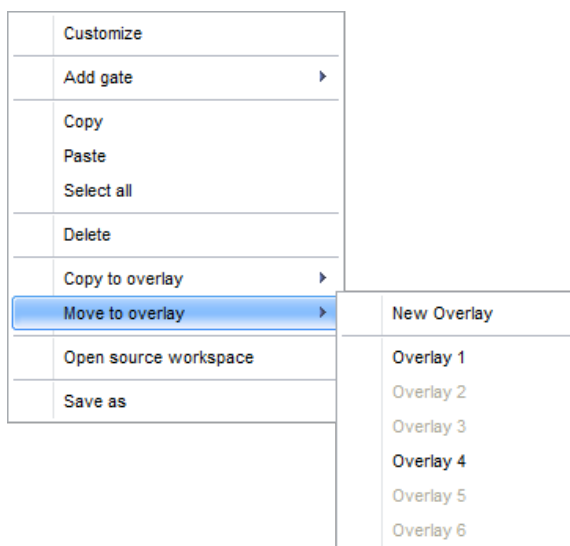
- ① Three plots selected as part of the multiselect are bounded by dark blue rectangles
- ② Plot that is not selected
- Alternatively, drag a selection rectangle using the mouse to select several plots within the same area.  
When drawing a selection rectangle, the selection process must be initiated outside a currently selected plot to avoid dragging the current plot to a new position.
- To deselect a single plot, click the selected plot as you hold down the **Ctrl** key. To deselect all plots, click an object other than a plot.

## Select all plots

- To select all plots in the **Overlays area**, press the **Ctrl+A** keys when an **Overlay plot** is selected.
- To select all **Gallery plots** for the selected **Overlays**, press the **Ctrl+A** keys when a **Gallery plot** is selected.

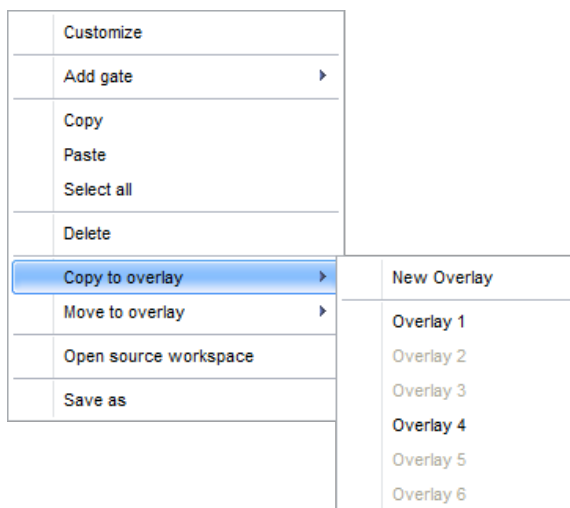
## Move and copy plots between Overlays

- To move a plot, right-click the **Gallery plot** to open the **Gallery plot context menu**, then select **Move to overlay** (“Gallery plot context menu options” on page 256).



Alternatively, plots can be moved from one **Overlay** to another by the cutting and pasting, using the standard **Ctrl+X** (cut) and **Ctrl+V** (paste) key combinations.

- You can copy a plot from one **Overlay** to another by using **Copy** and **Paste** options or the **Copy to overlay** option on the **Overlay plot context menu** or by using the standard **Ctrl+C** (copy) and **Ctrl+V** (paste) key combinations.





## Copy plot images

You can copy images of the selected plots in the **Overlay view** to the clipboard for pasting in other applications. This is accomplished by using the **Copy** option on the **Home ribbon tab** (“Home tab” on page 74) or using the standard **Ctrl+C** key combination.

## Resize plots

- **Overlay plots** in the **Overlay area** and **Gallery plots** in the **Gallery area** have a uniform size. You can resize **Overlay plots** and **Gallery plots** independently.
- Turn the **mouse scroll wheel** while holding down the **Ctrl** key to resize the **Overlay plot** or the **Gallery plots**, when dragging is not currently taking place.
- If an **Overlay plot** or a **Gallery plot** is selected, you can use the **Size Slider** in the lower right on the **Application Status Bar** (“Size slider” on page 70) to resize the selected plot.

## Overlay plot links to Source plot

### Link to live data of the Source plot

- **Overlay plots** are automatically updated to reflect the changes made to them in the **Workspace**, except for the **gates** created on **Source plots** in the **Workspace area**, which are not displayed on the **Overlay plots**.
- If a single-parameter plot is changed to a dual-parameter plot or vice versa, the **plot** is removed from its respective **Overlays**.

### Source plot not available

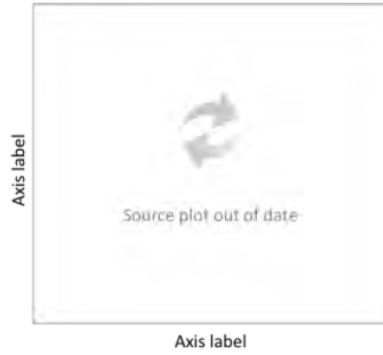
- When **Source plots** are deleted from the **Workspace** or a **Sample** is deleted, their corresponding plots are removed from the **Gallery** and the **Overlay area**.

### Empty Overlays

- If all **plots** in the same **Gallery** are deleted, the **Overlay** itself is automatically deleted.

## Auto-Refresh

- When **Auto-Refresh** is selected on the **Home ribbon tab** (“Home tab” on page 74), all plots and statistics in the **Overlays tab** are updated automatically in the background as the **Workspace** is modified.
- If **Auto-Refresh** is disabled or the system has not yet retrieved the image, the **Overlays** only show data from source locations that are up-to-date.
- When **Auto-Refresh** is disabled, the **Gallery plots** for the source locations that must be refreshed show the **out-of-date watermark** that states “**Source plot out of date**” on the plot.



The **Gallery plots** for the source locations that do not contain data are left blank.



- Only **Gallery plots** containing data (no watermark) are shown in the **Overlay plot**.

# Overlay plot context menu

## Overview

The **Overlay plot context menu** is displayed when an area within the plot boundary is right clicked.

The **Overlay plot context menu** contains commands for customizing the Overlay, adding gates to Overlay plots, for cutting, copying, pasting, and deleting selected plots, for sending selected plots to a selected Overlay, and saving all selected items as a new graphics file.

Customize
Add gate ▶
Copy
Paste
Select all
Delete
Copy to overlay ▶
Save as

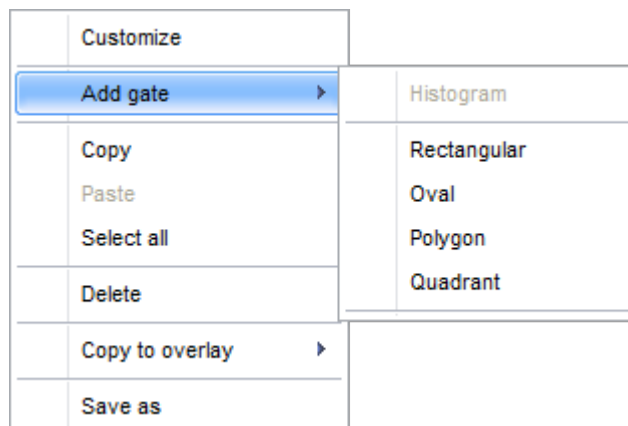
## Overlay plot context menu options

### Customize

- **Customize** opens the **Customize panel**, which enables you to modify the appearance of the selected **Overlay plot** as described in “Customize Overlay options” on page 474.
- If the panel is closed, selecting **Customize** opens the panel and brings it to the front of other open panels.

## Add gate

- **Add gate** creates a new gate in the default position on the currently selected plot. If multiple plots are selected, this option adds a gate to a single plot. The **Histogram** option is the only available for **Histogram plots**. **Rectangular**, **Oval**, **Polygon**, and **Quadrant** options are available only for **dual-parameter plots**.



- Selecting these options creates the selected gate type in the default position as described in “Create and delete gates” on page 145.
- The maximum number of gates allowed is 5, as described in “Gates” on page 240.
- The **Quadrant** option is unavailable if a **Quadrant gate** has been created on the current plot; all other options are unavailable after the gate limit has been reached.
- The gate set on the selected **Overlay plot** appears on all **Gallery plots** for the current **Overlay** and on the **Overlay plot**.

## Copy

**Copy** copies the selected plots to the clipboard, keeping the original plots in the **Overlays**.

## Paste

**Paste** inserts the objects currently in the clipboard onto the current **Overlay**.

## Select all

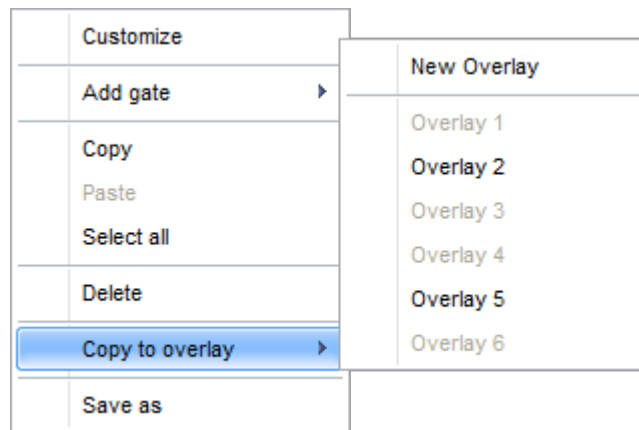
**Select all** selects all **Overlay plots**.

## Delete

**Delete** removes all selected **Overlay plots** and deletes all associated **Gallery plots**.

## Copy to Overlay

- **Copy to Overlay** sends a copy of the selected plot in the **Overlays area** to a selected **Overlay**, while retaining the original plot in the **Overlays area**.
- A new **Overlay** can be created by selecting **New Overlay** or by selecting an existing **Overlay** that contains plots with the same number of parameters.
- **Overlays** with non-matching number of parameters are listed in the **Overlay plot context menu**, but they cannot be selected.
- Similarly, if the selected number of plots causes an **Overlay** to exceed the allowable number of plots, that **Overlay** cannot be selected from the menu.



## Save as

- **Save as** opens the **File Save (Export) dialog** as described on “File Save (Export) dialog” on page 715, which enables you to save all selected plots as a new graphics file.
- The layout of the graphics file replicates the layout of the selected items on the current **Overlays view**, and the plot contents are based on the options selected in the **Export Options** tab of the **Options dialog** (see “Export Options” on page 662).
- The selection area for saving can only cover one page.
- The available image file formats are **PNG, BMP, GIF, JPG, TIF, and EMF** (Windows™ Metafile). The default format is **PNG**.

## Gallery plot context menu

### Overview

The **Gallery Plot context menu** is displayed when any area within the Gallery plot boundary is right clicked.

The **Gallery Plot context menu** contains the commands for customizing the Overlay, adding gates to Overlay plots, for cutting, copying, pasting, and deleting selected plots, for sending selected plots to a selected Overlay, and saving all selected items as a new graphics file.

Customize
Add gate ▶
Copy
Paste
Select all
Delete
Copy to overlay ▶
Move to overlay ▶
Open source workspace
Save as

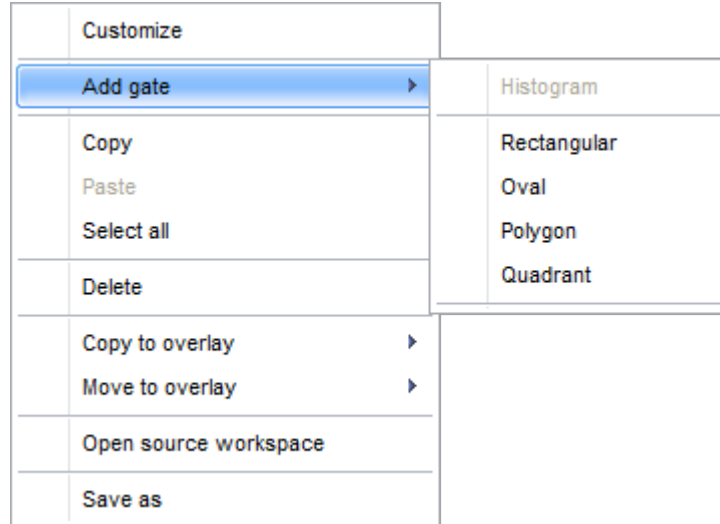
### Gallery plot context menu options

#### Customize

- **Customize** opens the **Customize panel**, which enables you to modify the appearance of the selected **Gallery plot** as described on “Customize Overlay options” on page 474.
- If the panel is closed, selecting **Customize** opens the panel and brings it to the front of other open panels.

#### Add gate

- **Add gate** creates a new gate in the default position in all gallery plots as described in “Create and delete gates” on page 145.
- The **Histogram** option is only available for **Histogram plots**.  
**Rectangular**, **Oval**, **Polygon**, and **Quadrant** options are available only for **dual-parameter plots**.
- The maximum number of gates allowed is 5, as described in “Gates” on page 240.
- The **Quadrant** option is unavailable if a **Quadrant gate** has been created on the current plot; all other options are unavailable after the gate limit has been reached.



## Copy

**Copy** copies the selected plots to the clipboard, keeping the original plots in the **Overlays**.

## Paste

**Paste** inserts the previously copied **Gallery plots** onto the current **Overlay plot**. The plots are appended at the end of the existing plots.

## Select all

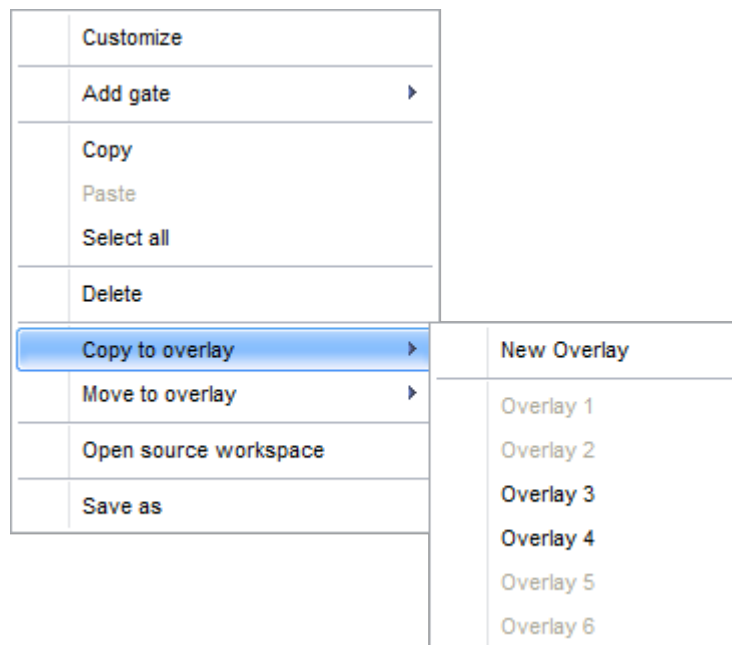
**Select all** selects all **Gallery plots** for the current **Overlay plot**.

## Delete

**Delete** removes selected **Gallery plots** and deletes them from the **Overlay**.

## Copy to Overlay

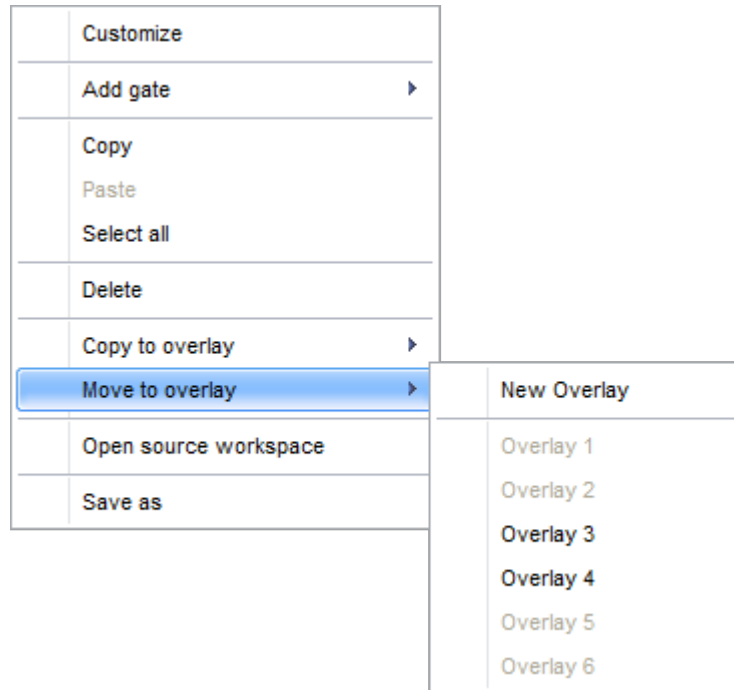
- **Copy to Overlay** sends a copy of selected plots in a **Gallery** to a selected **Overlay**, while retaining the original plots in the **Gallery**.
- A new **Overlay** can be created by selection of **New Overlay** or send the plots to an existing **Overlay** that contains plots with the same number of parameters.
- **Overlays** with non-matching number of parameters are listed in the **Gallery plot context menu**, but they cannot be selected.
- Similarly, if the selected number of plots causes an **Overlay** to exceed the allowable number of plots, that **Overlay** cannot be selected from the menu.



## Move to overlay

- **Move to overlay** moves the selected plots within a **Gallery** to a selected **Overlay**. The **Gallery** does not retain the plot that is moved.
- A plot can be moved to either a **New Overlay** or an existing overlay.
- **Overlays** with non-matching number of parameters are listed in the **Gallery plot context menu**, but they cannot be selected.
- Similarly, if the selected number of plots causes an **Overlay** to exceed the allowable number of plots, that **Overlay** cannot be selected from the menu.





## Save as

- **Save as** opens the **File Save (Export) dialog** as described on “File Save (Export) dialog” on page 715, which enables you to save all selected plots as a new graphics file.
- The layout of the graphics file replicates the layout of the selected items on the current **Overlays view**, and the plot contents are based on the options selected in the **Export Options** tab of the **Options dialog** (see “Export Options” on page 662).
- The selection area for saving can only cover one page.
- The available image file formats are **PNG**, **BMP**, **GIF**, **JPG**, **TIF**, and **EMF** (Windows™ Metafile). The default format is **PNG**.

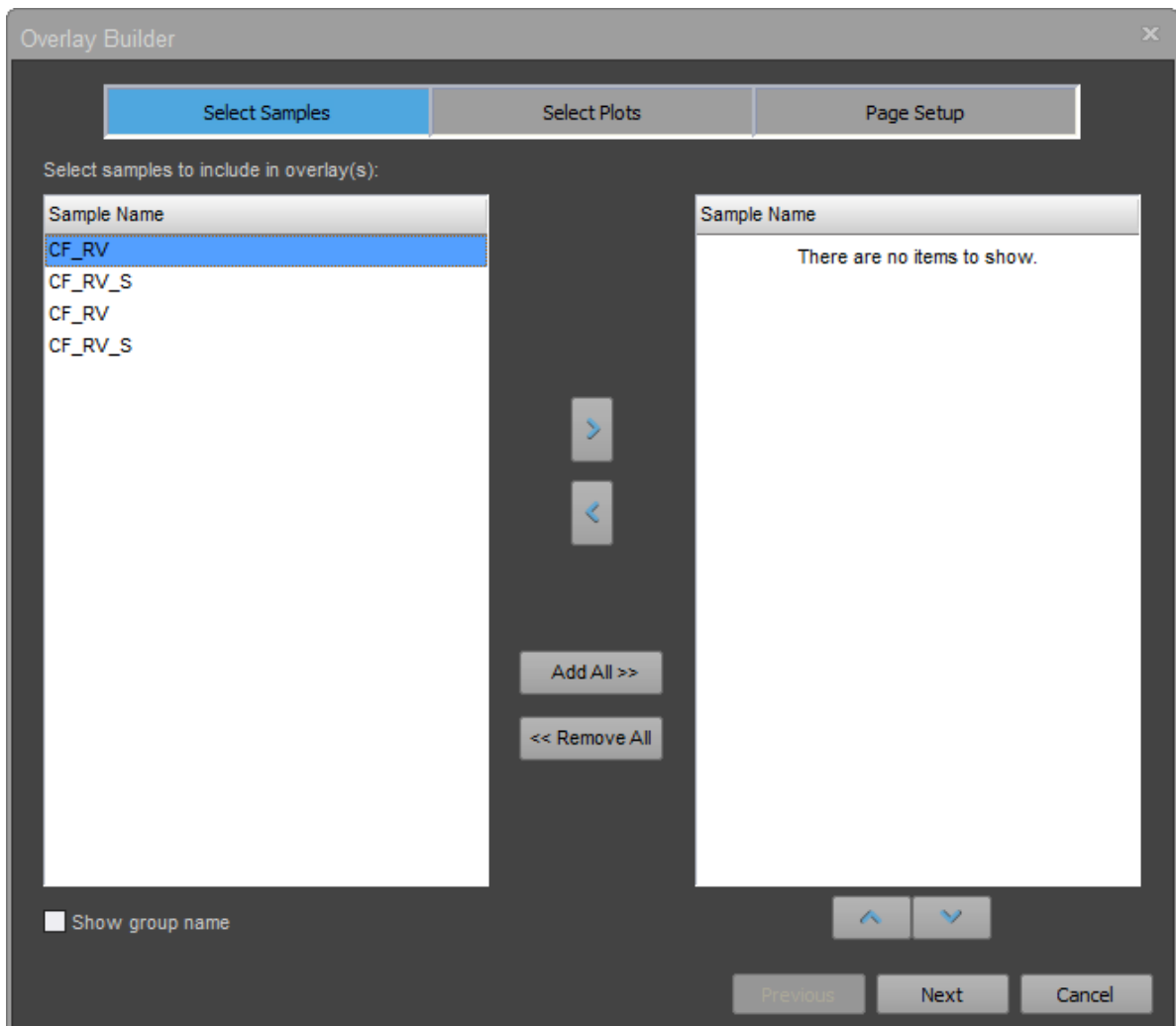
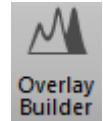
# Overlay Builder

## Overview

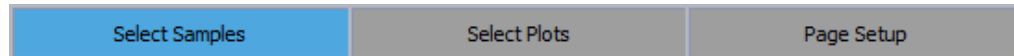
The **Overlay Builder** tool guides you through the process of creating new **Overlay plots** in a single **Experiment**. The **Overlay Builder** enables you to select samples and plots for a new **Overlay** and determine how to display the selected plots.

**Note:** Overlay of plots from different Experiments is not permitted.

- To open the **Overlay Builder**, click the **Overlay Builder** button on the **Overlay ribbon tab** (“Overlay tab” on page 110). **Overlay Builder** is only enabled when at least one plot exists on the Experiment Workspace and the instrument is not acquiring.



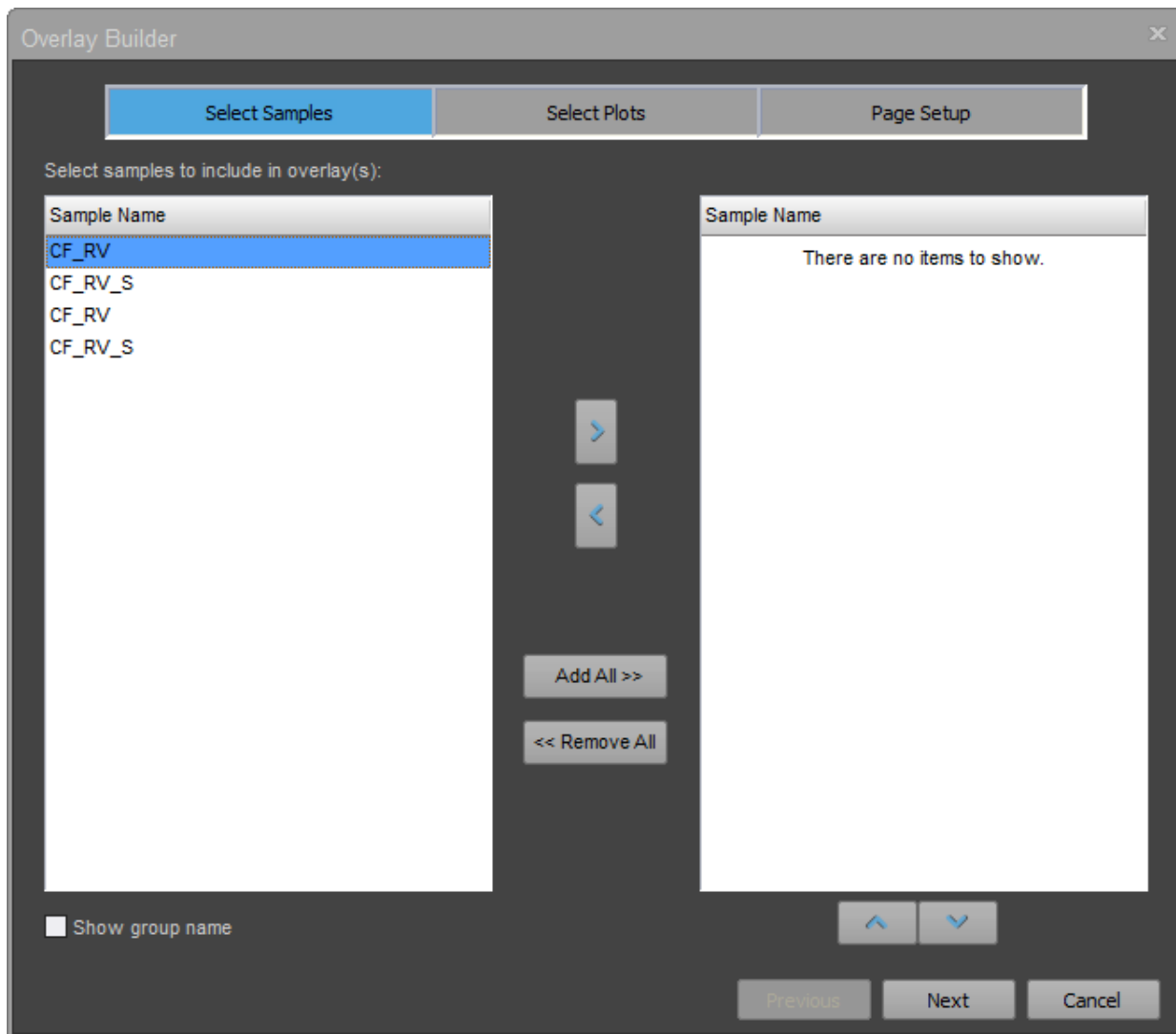
- The **Overlay Builder** consists of three tabs, listed below, that are displayed sequentially to guide you through the process of building a new Overlay.
  - **Select Samples** (“Select and order samples” on page 262)
  - **Select Plots** (“Select Plots” on page 264)
  - **Page Setup** (“Page Setup” on page 266)
- The top portion of the **Overlay Builder** contains the tabs, where the current tab is highlighted.



- **Next** moves to the next tab of the **Overlay Builder**.  
The **Next** button is not displayed on the last tab of the builder (**Page Setup** tab), where it is replaced by the **Finish** button.
- **Previous** returns to the previous tab of the builder.  
The **Previous** button is not active on the first tab of the builder (**Select Samples** tab).
- **Finish** closes the **Overlay Builder** and creates a new **Overlay** based on the selections made in the **Overlay Builder**.
- **Cancel** closes the **Overlay Builder** without creating **Overlay plots**. Any changes made with the builder are discarded.  
The **Cancel** button is available on all pages of the builder.
- The default title for a new Overlay is **Overlay n**, where **n** is a numerical suffix (first available integer) that ensures that the title is unique.  
**Gallery plot titles** show the **Sample name** by default.

## Select and order samples

The **Select Samples** tab of the **Overlay Builder** enables you to select samples for inclusion in the New Overlay and reorder the selected samples, if desired.



## Sample lists

- The **Available Samples** panel on the left contains a list of all **Samples** in the current **Experiment** based on the Experiment hierarchy.  
By default, this list only displays the **Sample name**.  
Select the **Show group name** option to show the **Group name** in front of the **Sample name** (e.g. group1\sample1).
- The **Selected Samples** panel on the right contains a list of the **Samples** that are selected to be displayed in the **Overlay plots**. By default, this list is empty.
- **Sample names** only appear in one list at a time.
- Use the **arrow** buttons (“Arrow buttons” on page 263) located between the two panels to transfer samples between the two lists.
- You can select multiple **Samples** in a list simultaneously using one of the following methods:
  - To select non-consecutive **Samples** in a list, press and hold down the **Ctrl** key, then click each **Sample** that you want to select.
  - To select a consecutive group of **Samples** in a list, click the first **Sample**, press and hold down the **Shift** key, then click the last **Sample**.
  - Use the **Ctrl+A** key combination to select all **Samples** in the currently active list.
- Vertical and horizontal scroll bars appear in each list if needed.

## Arrow buttons

- Click the **right arrow** to transfer all selected samples from the **Available Samples** panel to the **Selected Samples** panel.  
The **right arrow** is only enabled when one or more **Samples** are selected in the **Available Samples** list.
- Click the **left arrow** to transfer all selected **Samples** from the **Selected Samples** panel to the **Available Samples** panel.  
The **left arrow** is only enabled when one or more **Samples** are selected in the **Selected Samples** list.

## Add All

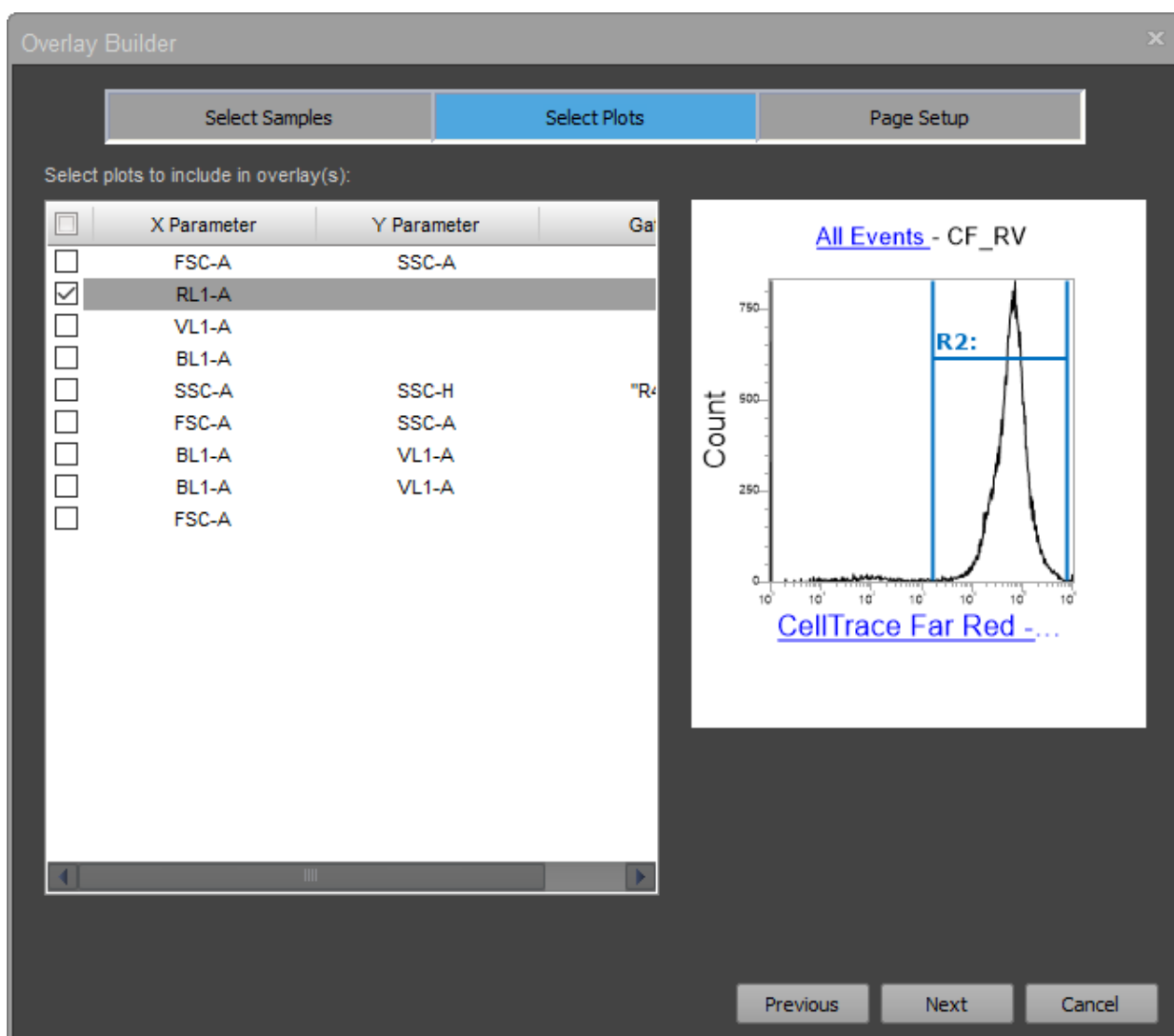
- Click **Add All** to transfer all **Samples** from the **Available Samples** panel to the **Selected Samples** panel.

## Remove All

- Click **Remove All** to transfer all **Samples** from the **Selected Samples** list to the **Available Samples** panel.

## Select Plots

The **Select Plots** tab enables you to select plots to display them on an Overlay plot.



- The **Select Plots** tab contains the **Plot Selection Table** (left panel) and a **preview image** of the currently selected plot (right panel).

## Plot selection table

- The **Plot Selection Table** lists all plots present on the Experiment-level Workspace of the current Experiment in the order they were created.

<input type="checkbox"/>	X Parameter	Y Parameter	Gate
<input type="checkbox"/>	FSC-A	SSC-A	
<input checked="" type="checkbox"/>	RL1-A		
<input type="checkbox"/>	VL1-A		
<input type="checkbox"/>	BL1-A		
<input type="checkbox"/>	SSC-A	SSC-H	"R4"
<input type="checkbox"/>	FSC-A	SSC-A	
<input type="checkbox"/>	BL1-A	VL1-A	
<input type="checkbox"/>	BL1-A	VL1-A	
<input type="checkbox"/>	FSC-A		

The **Plot Selection Table** has columns for the **X Parameter**, **Y Parameter**, and parent **Gate** name columns, and an extra **Select** column to the left of the plot title used for selecting the plots of interest.

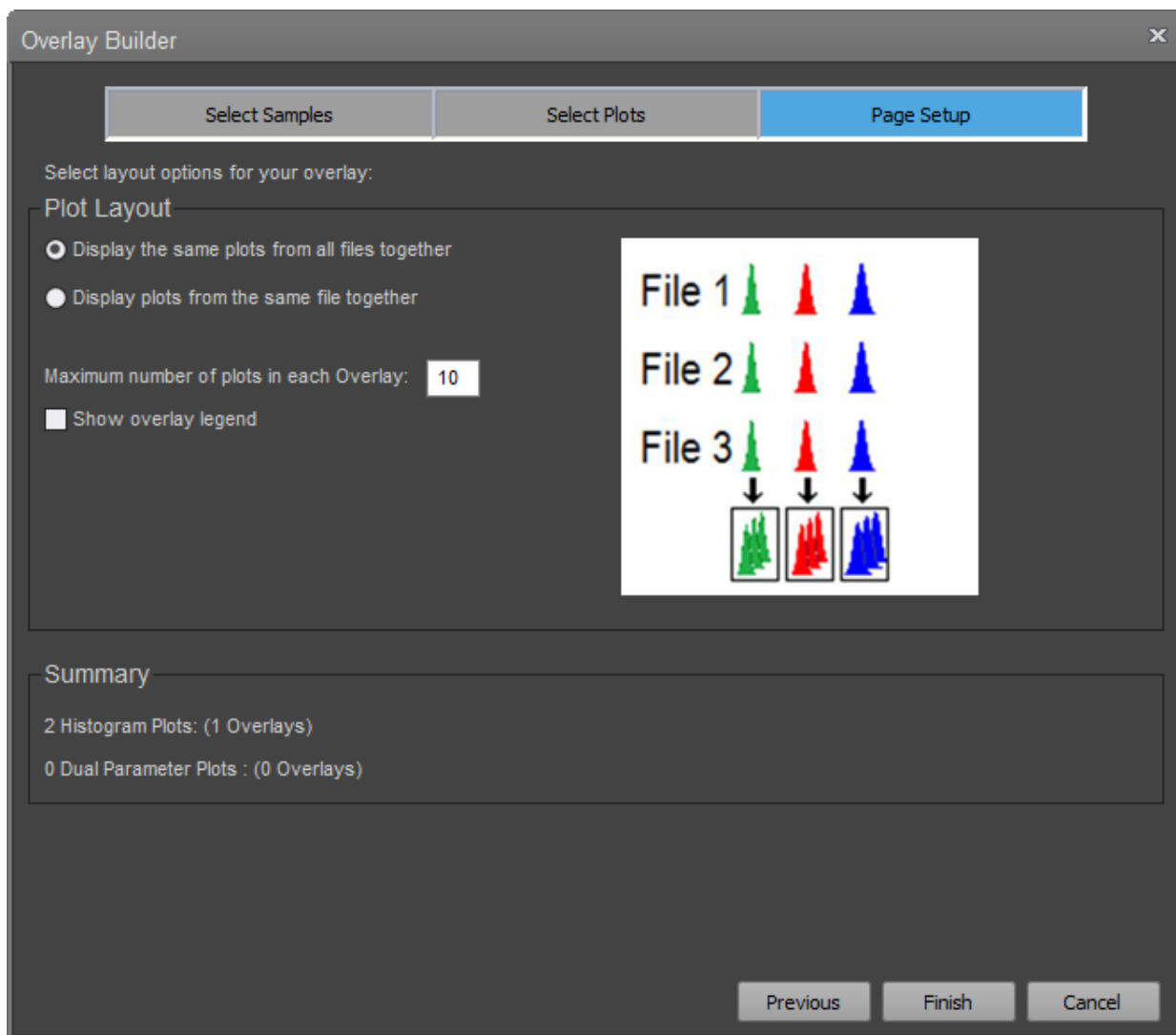
- To add one or more plots to an Overlay plot, select the relevant **checkboxes** next to the plot of interest.
- To add all plots listed in the plot selection list, click the **select column heading checkbox**.
- Deselect the **select column heading checkbox** to deselect all individual plots.
- If one or more of the plot checkboxes are deselected, the select column heading checkbox is also unchecked.
- By default, no plots are selected for overlay and all checkboxes are deselected.
- Vertical and horizontal scroll bars appear in the list if needed.

## Plot preview

- The **Plot preview** section of the **Overlay Builder** uses plots from the Experiment workspace regardless of what workspace is active.
- Click a single row in the plot selection list to highlight the selected row and show a preview of the selected plot in the preview area.
- The preview shows the selected plot with data from the first selected sample together with any gates and labels that are present on the plot.
- You cannot format plots or modify gate positions on the preview image.

## Page Setup

**Page Setup** enables you to set the display options for the selected plots.

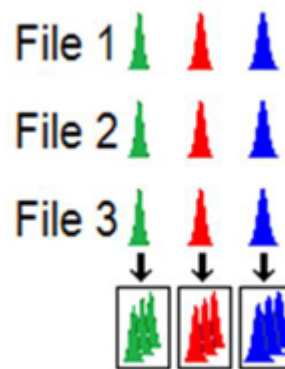


- The **Plot Layout** section contains options to control how selected plots from selected files are added to new Overlay plots.

### Display the same plots from all files together

- The **Display the same plots from all files together option** is the default option. It creates **Overlay plots** that include data from all samples (multiple files) in an Experiment.

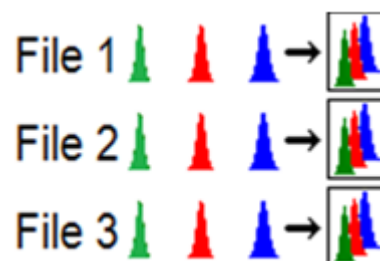




- When the **Display the same plots from all files together** is selected, **Overlay plots** are created according to the following rules:
  - Selected **Histogram plots** are placed in an **Overlay plot** based on the **X-axis parameter**. Each parameter is contained in a different **Overlay plot**.  
If only one **Histogram plot** is selected for **Overlay**, only one **Histogram Overlay plot** is generated.  
If multiple plots are selected, multiple **Histogram Overlay plots** are generated.
  - If the number of plots set in the **Maximum number of plots in each Overlay** (see “Maximum number of plots in each Overlay” on page 268) is exceeded at any time, extra **Overlay plots** are created to contain the remaining single-parameter **Histogram plots** from each file.
  - All selected **dual-parameter plots** are contained in an **Overlay plot** based on like **X-axis** and **Y-axis parameters**. Each parameter combination is placed in a different **Overlay plot**.  
If only one dual parameter plot is selected for **Overlay**, only one **Overlay plot** of dual parameter data is generated.  
If multiple plots are selected, multiple dual parameter **Overlay plots** are generated.
  - If the number of plots set in the **Maximum number of plots in each Overlay** is exceeded at any time, extra **Overlay plots** are created to contain the remaining dual-parameter **Histogram plots** from each file.

### Display plots from the same file together

- When the **Display plots from the same file together** option selected, each **Overlay plot** generated contains plots from a single Sample (one FCS file) in the Experiment and not multiple Samples. This option is useful if fluorescence emission is compared between parameters, for example spillover of FITC into PE.



- When **Display plots from the same file together** is selected, **Overlay plots** are created according to the following rules:
  - All single-parameter **Histograms** specific for the Sample are placed in a single **Overlay plot**.
  - If the number of plots set in the **Maximum number of plots in each Overlay** (see “Maximum number of plots in each Overlay” on page 268) is exceeded at any time, extra **Overlay plots** are created to contain the remaining single-parameter **Histogram plots** from the same file.
  - All selected **dual-parameter plots** specific for the Sample are placed in a single **Overlay plot**.
  - If the number of plots set in the **Maximum number of plots in each Overlay** is exceeded at any time, more **Overlay plots** are created to contain the remaining dual-parameter **Histogram plots** from the same file.

### Maximum number of plots in each Overlay

- **Maximum number of plots in each Overlay** lets you specify the maximum number of plots to be added to an **Overlay plot** by the **Overlay Builder**. If the requested plots exceed this number, new **Overlay plots** are created as needed.
- The default number of plots in an Overlay is 10 plots.  
The minimum number of plots is 1.  
The maximum number of plots is 400.  
If you attempt to enter a number outside this range, the number is automatically adjusted to the closest limit.
- Only numeric characters are allowed.
- If the number entered is greater than the number of samples or plots, then the Overlay is created with the number of samples or plots.

### Show overlay legend

- If the **Show overlay legend option** is checked, **Overlay plots** are created with the legend panel displayed as described in “Legend” on page 238.

### Summary

- **Summary** shows the number of **Source plots** and **Overlay plots** created for both **Histogram** and **dual parameter plots**.

## Print overlays

The print command (**Ctrl+P**) and previews are enabled when either the **Overlay plot** or **Gallery view** is in focus.

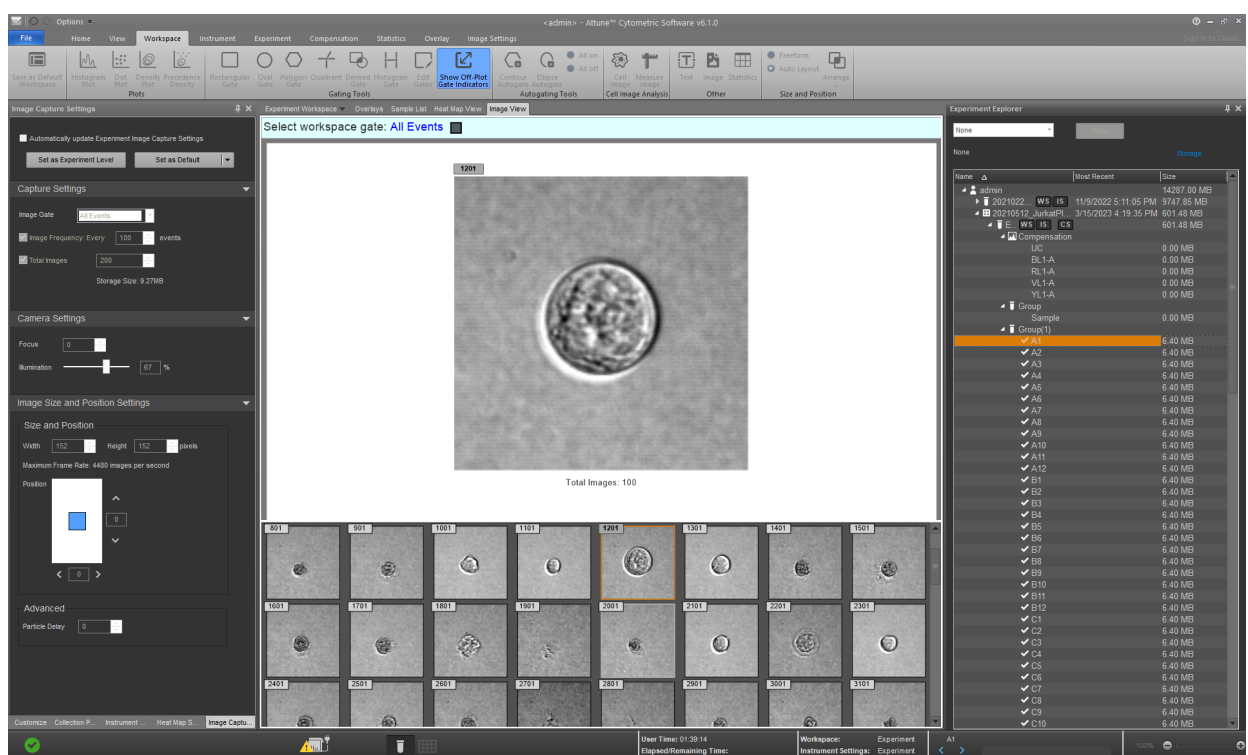
- The **Overlays** are printed in the order they appear on the **Overlays tab** using the current page layout settings (orientation and paper size).
- **Gallery plots** are not printed.
- The zoom setting of the **Overlays** affects the printed size of the overlay plot. The number of **Overlays** per page is based on the zoom and the page layout settings.
- Before the **Overlays** are printed, the **Statistics** are refreshed.

## Overview

**Image View** enables you to view the cell images that are collected from a Sample during acquisition. In addition, you can:

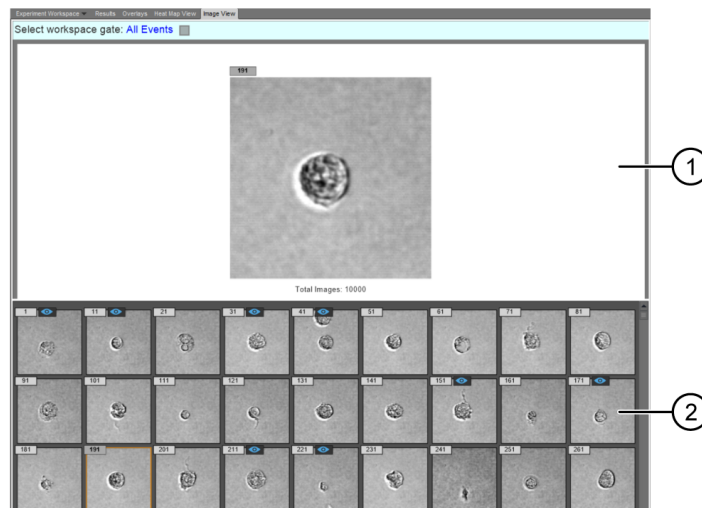
- Select from which population the images are captured using an **Image gate** (“Image capture gate” on page 276)
- Backgate displayed images to identify the associated events on the corresponding plot in Workspace view (“Image backgating” on page 278)
- Measure the area of imaged cells (“Measure image tool” on page 280)

**Note:** **Image view** is only available when using an Attune™ CytPix™ Flow Cytometer or if the current Experiment was created with an Attune™ CytPix™ Flow Cytometer.



## Image view tab

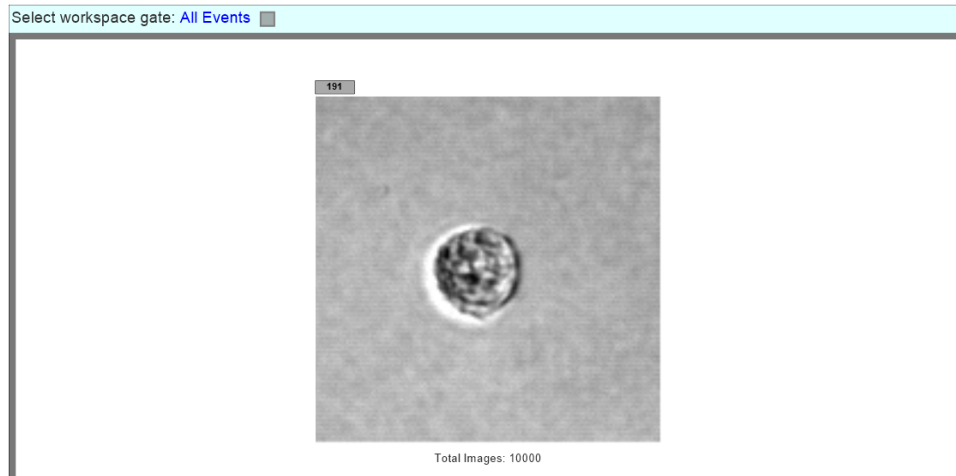
- **Image view** tab is displayed in the Application area on the **Main Application Workspace** (“Main application workspace” on page 56).
- The **Image view** tab is only visible when the **Attune™ desktop** is in view. It is not visible on the **Main menu**, **Login screen**, **Performance test views**, or **Filter configuration views**.
- By default, the **Image view** tab is docked to the center application window.  
The **Image view** tab can be undocked, center docked (with main application view windows), left docked, or right docked. The tab’s docked state, size, and position persists as part of the user settings.
- You can toggle the **Image view** on and off using the **View Images** toggle button in the **Image Settings** ribbon tab (“Image Settings tab” on page 112).
- **Image view** tab consists of an **Active image area** and an **Image gallery** that contains all the captured images.
- The **Active image area** is separated from the **Image gallery** using a splitter window control that enables resizing of the **Active image** and the **Image gallery** areas.



- ① Active image
- ② Image gallery

## Active image

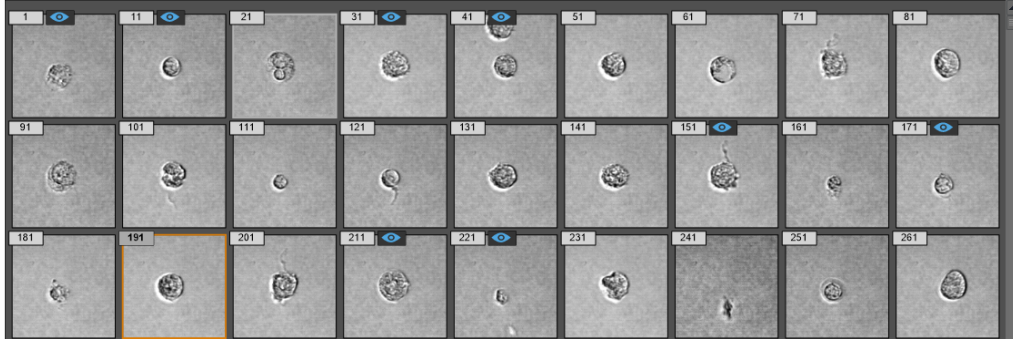
The **active image** is displayed in the center of the main active image area.



- By default, the active image area displays the first image that is acquired. If the active sample or the image gate does not contain any images, the image container displays the message “No images are available”.
- The active image maintains its aspect ratio and fills the available size of active image area when the active image is resized.
- The active image allows additional controls using the **Customize** panel, right-click **context menu options**, and **image analysis tools**.
- The active image area in **Image View** contains a **hyperlink** that enables you to select a gate from the Workspace to filter the images displayed in the **Image View** tab (see “Image capture gate” on page 276).

## Image gallery

The **Image gallery** displays all the images that are contained in a selected gate for the active workspace.



- The gallery displays up to 2000 images, which are arranged in the image acquisition event order. When the splitter window is adjusted to change the area available to the **Active Image** and the **Image Gallery**, the number of images per row in the **Image Gallery** is dynamically updated based on the available area. The minimum number of images per row is 1.
- Only images that are in view are loaded into memory. While images are being loaded, a refresh image is displayed in place of the captured image. After the captured image has been read from the disk, the image replaces the refresh image in the gallery.



- The gallery area contains vertical scroll bars to allow more images to be visible in rows.
- The images shown in the gallery show the **event ID** in the top left of the image.
- In the event an image is missing, the **parameter not available** icon serves as a placeholder and the **event ID** is displayed in the top left of the image.



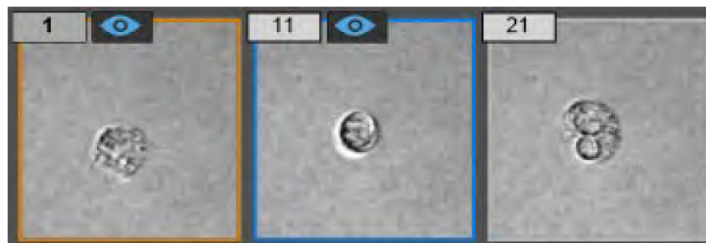
## Navigation and selection

- You can navigate the **Image gallery** using the **up**, **down**, **left**, and **right arrow** keys on the computer keyboard.
- **Page up** and **page down** keys scroll the view by one page up or down.
- The mouse scroll wheel moves the view up and down as a line per click of the mouse.
- You can select any image in the gallery to make that image the **active image**.

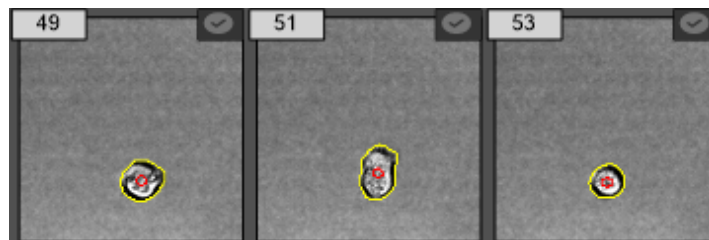
The **active selected image** is indicated with an orange border (image 1 in the example below).

A **selected image that is not the active image** is indicated with a blue border (image 11 in the example).

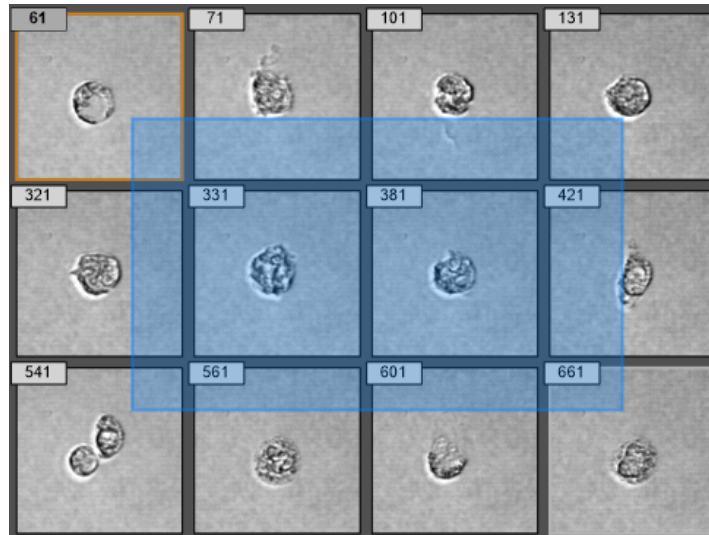
**Images that have been backgated** (“Image backgating” on page 278) are indicated with an **eye** icon (images 1 and 11 in the example).



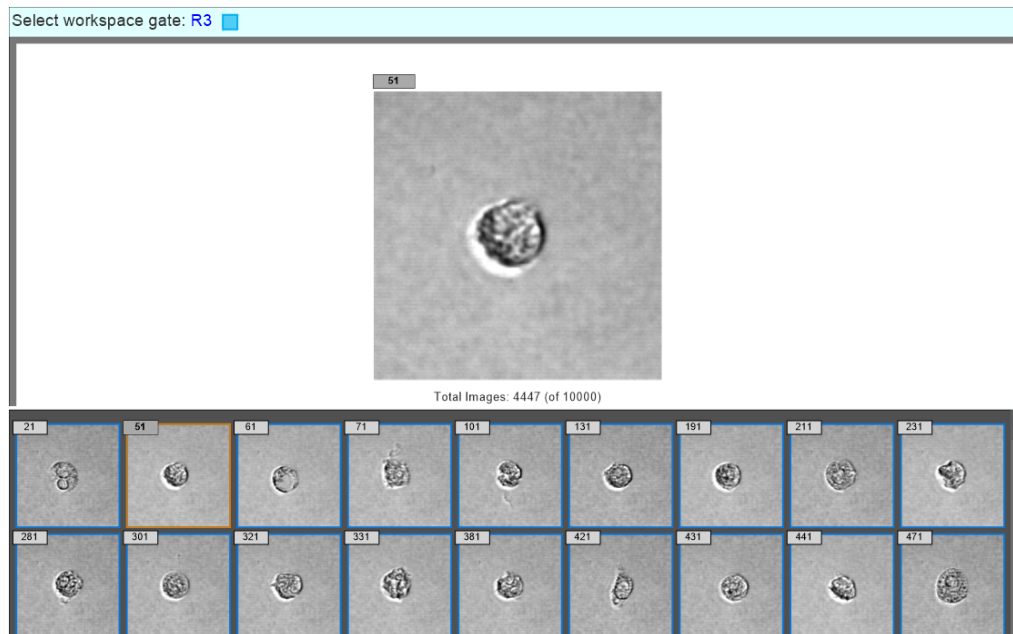
- **Images that have been processed** are indicated with a check icon (images 49, 51, and 53 in the example below).



- You can select multiple images using one of the following methods:
  - Left-click and drag to select all images within the defined rectangle.

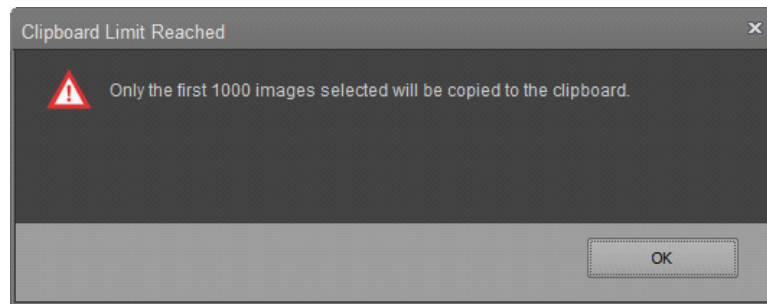


- Click as you hold **Ctrl** to select or deselect non-contiguous images.
  - Press **Ctrl+A** to select all images in the gallery.
- When multiple images are selected, the images that are selected are indicated with a blue border. The **active image** in the selection is still indicated with an orange border (image 51 in the example below).



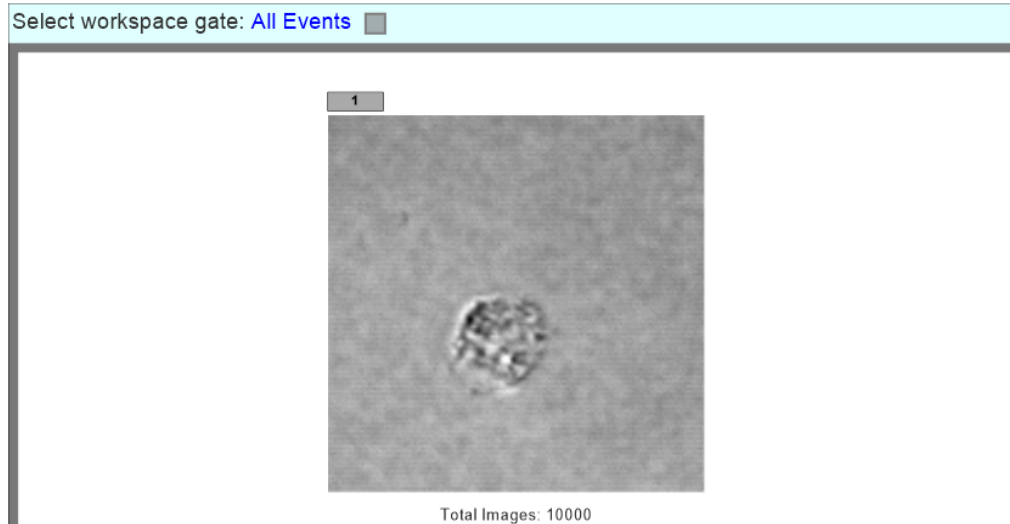


- Only 1000 images can be selected to the clipboard or sent to the Workspace. If more than 1000 images are selected, the software displays the **Clipboard Limit Reached** dialog, which states that only the first 1000 images selected will be copied to the clipboard.

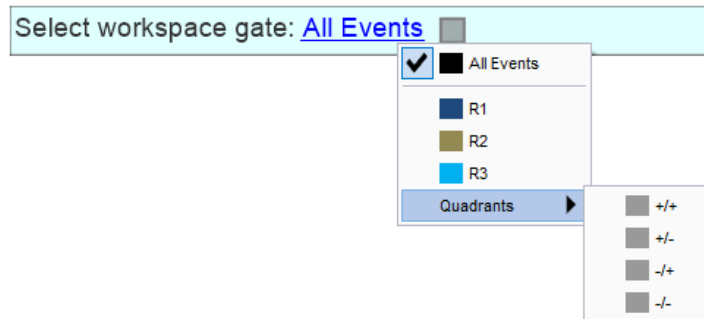


## Image capture gate

The **active image area** in the **Image view** contains a **Select workspace gate** hyperlink that enables you to select from which population the images are captured (**Active image** and **Image Gallery images**) using an **Image capture gate**.

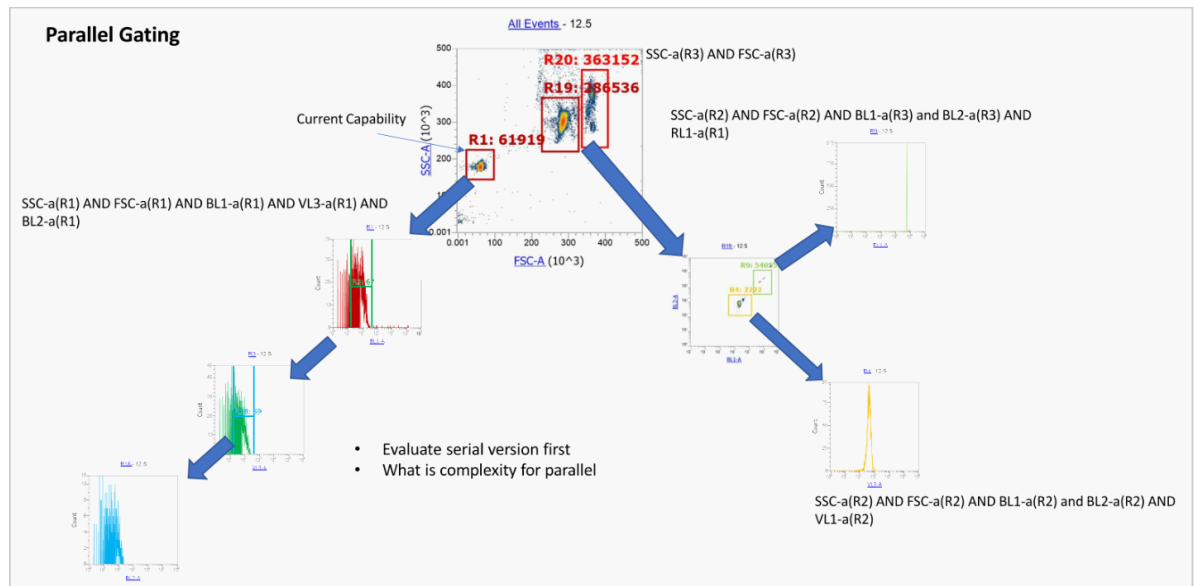


- When the **Select workspace gate** hyperlink is selected, the **dynamic gate context menu** is displayed.

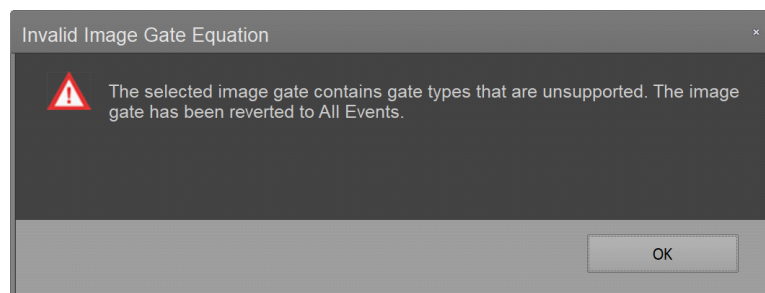


- The **gate context menu** displays a list of all available gates from the active Workspace and includes the **All Events** option as well as any derived gates.
- The **gate context menu** only shows gates that are valid.
  - **Image capture gate types** are limited to **rectangle**, **histogram**, and **non-bent quadrant gates**.
  - **Polygon**, **bent quadrant**, and **elliptical gates** are **not** supported.
  - **Derived/Logic gates** are supported as long as gate terms do not contain unsupported gate types.

- The **image capture gate** function supports up to 8 parallel gating hierarchies (i.e., **ORed gates**) with up to 20 levels.



- When you select a gate, only the images of events that are contained in the selected gate are displayed in the **Image view**.
- Compensation** is **not** applied to the data at the point where image decision is made.
- If the source gate is moved, the images in the image filter are updated to reflect the event images contained in the moved gate.
- If the source gate is deleted, the image filter reverts to **All Events**.
- If any gates in the gating hierarchy are changed to an invalid gate type (i.e., **polygon**, **autogate**, etc.), the **image capture gate** reverts to **All Events**. In this case, the software displays the **Invalid Image Gate Equation** dialog, which states that an invalid gate type was used and that the image capture gate will revert to **All Events**.



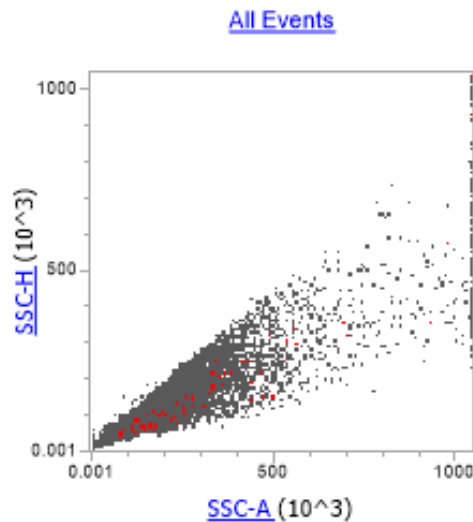
- The **image capture gate** function is available only if the selected sample's listmode specification contains both the **event** and **imageflag** parameters. Otherwise, it reverts to **All Events**.
- When available, the **image capture gate** function applies to both a recorded file and empty sample as long as the sample is contained in an Experiment that supports imaging and the active instrument model is set to an imaging cytometer model (i.e., an Attune™ CytPix™ Flow Cytometer).

## Image backgating

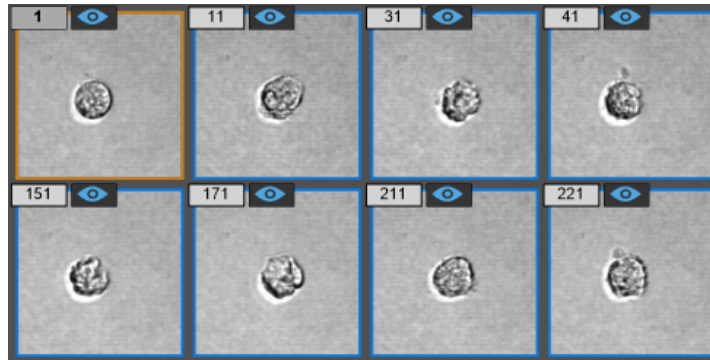
**Image backgating** enables you to identify the event that is associated with the captured cell image on a **Dot plot** or a **Precedence Density plot** in Workspace view (“Image backgating” on page 179).

**Note:** **Image backgating** is only available when using an Attune™ CytPix™ Flow Cytometer or if the active Experiment was created with an Attune™ CytPix™ Flow Cytometer.

- To backgate images in **Image View**, right-click the **cell image** you want to backgate in **Image View**, then select **Show Image on Plots**. You can select the **Active image** or any image from the **Image Gallery**.  
To backgate multiple images, select the images you want to backgate in the **Image Gallery**, then right-click and select **Show Images on Plots**.
- By default, backgated events appear on the associated Workspace plot in red, but you can select a different color for the backgated events using the **Image Options** dialog (“Image Options” on page 673).



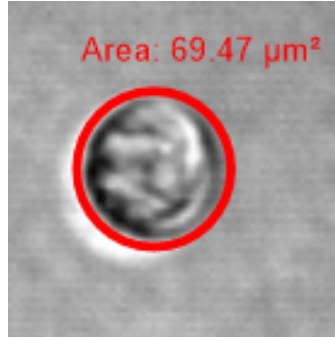
- The images that have an image backgate applied display an **eye** icon.



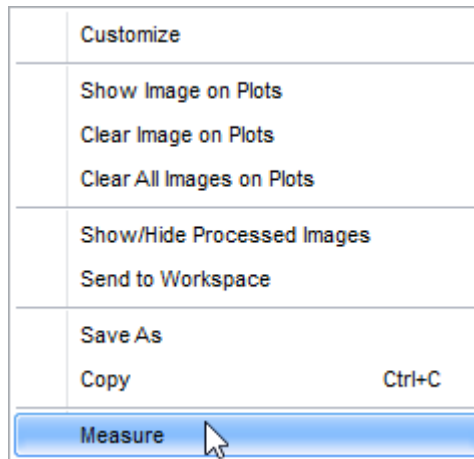
- To undo the backgating of an image, right-click the **backgated image**, then select **Clear Image on Plots**.  
To undo all backgates on a plot, right-click the **Active image** or on any image in the **Image Gallery**, then select **Clear All Images on Plots**.

## Measure image tool

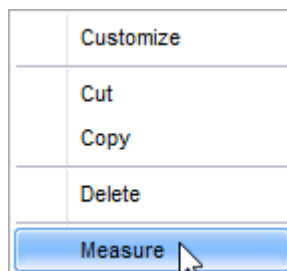
The **Measure image** tool enables you to measure the area of an imaged cell by drawing an ellipse on the active image that encircles the imaged cell.



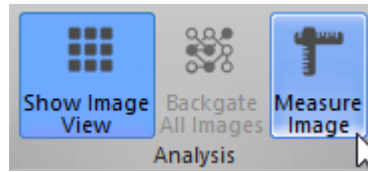
- You can select the **Measure image** tool by one these methods:
  - Right-click the **active image** in the **Image View** tab, then select **Measure** from the **active image context menu**.



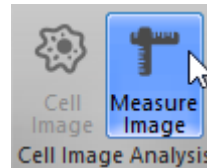
- Right-click a **cell image container** in the Experiment Workspace, then select **Measure** from the **cell image container context menu**.



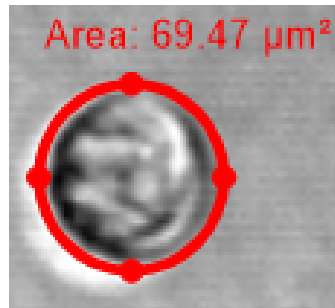
- Click **Measure Image** in the **Analysis** group of the **Image Settings** ribbon tab.



- Click **Measure Image** in the **Cell Image Analysis** group of the **Workspace** ribbon tab.



- You can resize the ellipse using the handles that become visible when the ellipse is selected to stretch and rotate it.
- To delete the measurement tool, click the ellipse, then select from the keyboard or the context menu.
- The measured area is based on a conversion factor of micrometers per pixel.



## Image view context menus

### Active image context menu

The **Active image context menu** is displayed when you right-click the **active image** in the **Image View** tab.

The **Active image context menu** contains these options:

Customize	
Show Image on Plots	
Clear Image on Plots	
Clear All Images on Plots	
Show/Hide Processed Images	
Send to Workspace	
Save As	
Copy	Ctrl+C
Measure	

- **Customize:** Opens the **Customize Image** panel, which enables you to adjust the **Brightness** and **Contrast** of the image, set **Scaling** options, and adjust **Mask Settings** (“Customize Image options” on page 484).
- **Show Image on Plots:** Identifies the event on the Workspace plots that is associated with the active image. By default, the backgated imaged event is shown in red.
- **Clear Image on Plots:** Removes the backgate associated with the image from the Workspace plots.
- **Clear All Images on Plots:** Removes all backgates associated with all images from the Workspace plots.
- **Show/Hide Processed Images:** Toggles between showing and hiding the **check** icon that identifies processed images in the **Image Gallery**.
- **Send to Workspace:** Sends the selected image to the active Workspace.
- **Save As:** Opens the **File Save (Export)** dialog, which enables you to save the selected image in the desired location (“File Save (Export) dialog” on page 715). The **Save As** option is disabled if acquisition is in progress.
- **Copy:** Copies the selected image to the windows clipboard. You can also press **Ctrl+C** on the keyboard to copy the selected image.
- **Measure:** Enables you to use the **Measure Image** tool with the selected image (“Measure image tool” on page 280).



## Image Gallery context menu

The **Image Gallery context menu** is displayed when you right-click the images displayed in the **Image Gallery** in the **Image View** tab.

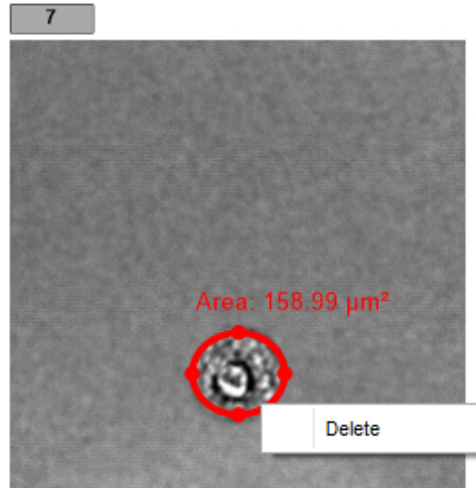
Customize
Show Selected Image(s) on Plots
Clear Selected Image(s) on Plots
Clear All Images on Plots
Show/Hide Processed Images
Send to Workspace
Save As
Copy <span style="float: right;">Ctrl+C</span>
Delete

The **Image Gallery context menu** contains the following options:

- **Customize:** Opens the **Customize Image** panel, which enables you to adjust the **Brightness** and **Contrast** of the image, set **Scaling** options, and adjust **Mask Settings** (“Customize Image options” on page 484).
- **Show Selected Image(s) on Plots:** Identifies the events on the Workspace plot that is associated with the selected images. By default, the backgated imaged event is shown in red.
- **Clear Selected Image(s) on Plots:** Removes the backgate associated with the selected images from the Workspace plots.
- **Clear All Images on Plots:** Removes all backgates associated with all images from the Workspace plots.
- **Show/Hide Processed Images:** Toggles between showing and hiding the **check** icon that identifies processed images in the **Image Gallery**.
- **Send to Workspace:** Sends the selected images to the active Workspace.
- **Save As:** Opens the **File Save (Export)** dialog, which enables you to save the selected images in the desired location (“File Save (Export) dialog” on page 715). The **Save As** option is disabled if acquisition is in progress.
- **Copy:** Copies the selected images to the windows clipboard. You can also press **Ctrl+C** on the keyboard to copy the selected images.
- **Delete:** Opens the **Confirm Deletion** dialog, which enables you to delete the selected images. The **Delete** option is disabled if acquisition is in progress.

## Measure image tool context menu

The **Measure image tool context menu** is displayed when you right-click the **image measurement tool**.



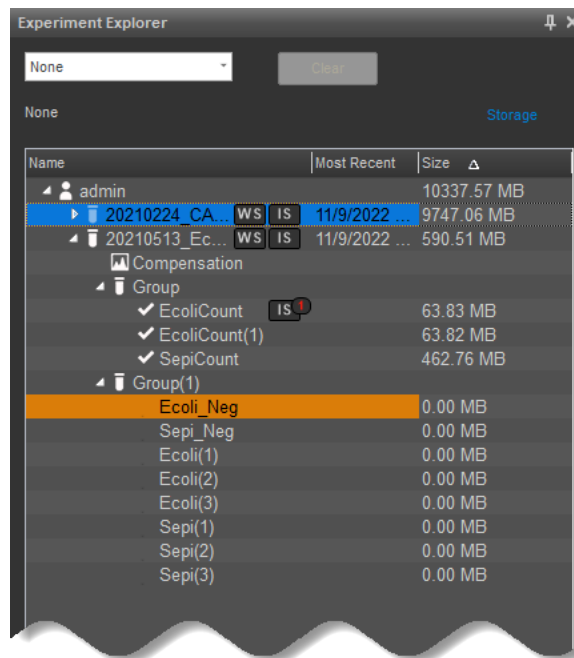
The **Measure image tool context menu** contains only the **Delete** command, which deletes the selected measurement tool.



# Experiment Explorer

## Overview

**Experiment Explorer** functions as an interface for creating, viewing, and managing Experiments.



**Experiment Explorer** is displayed both on the **Main Menu** (“Main Menu” on page 46) and the **Attune™ Desktop** (Chapter 3, “Attune™ Desktop”). By default, it is docked on the right side of the application window.

- **Files view** section (“Files view” on page 287) displays the user folders, which contain all **Experiment** files in a hierarchal view.
- The **Files view** section can contain multiple folders, each with multiple **Experiments**.
- Each **Experiment** can have more than one **Group**, each with one or more **Samples**.
- Each **Sample** is associated with its own FCS file.








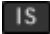



---

**Note:** When the Attune™ Cytometric Software is in the **SAE** mode, you can show application **object IDs** in the **Experiment Explorer** (see “Display ID” on page 299).

---

## Experiment Explorer icons

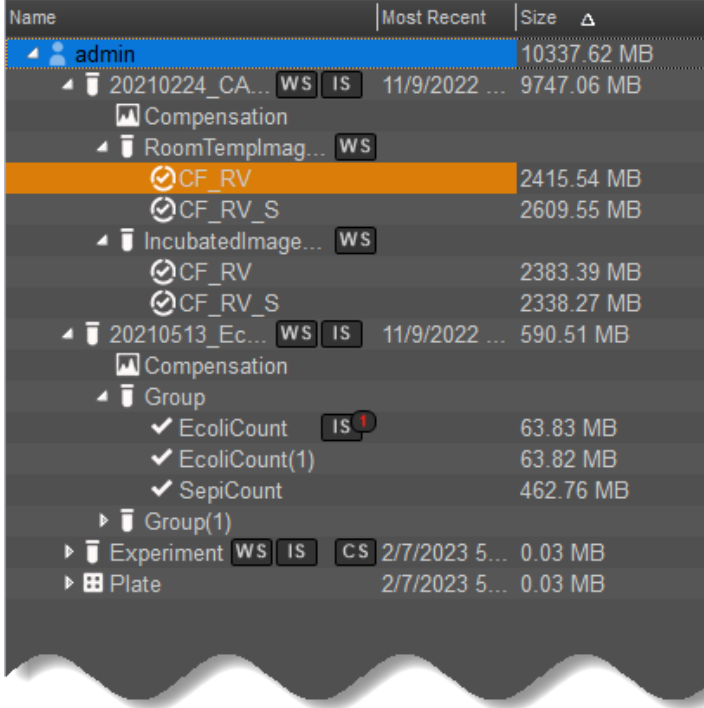
Experiment Explorer uses the following icons to indicate the various elements and their conditions in the Files view.

Icon	Indication	Icon	Indication
	User		Compensation
	Plate Experiment node		Read-only compensation
	Tube Experiment node		Workspace badge
	Group node		Instrument Settings badge
	Sample with FCS data		Compensation Settings badge
	Sample with FCS and Image Processing data		

# Files view

## Organization

Files view shows all Experiments and Samples generated by the current user or the Samples selected as a result of a filter being applied.

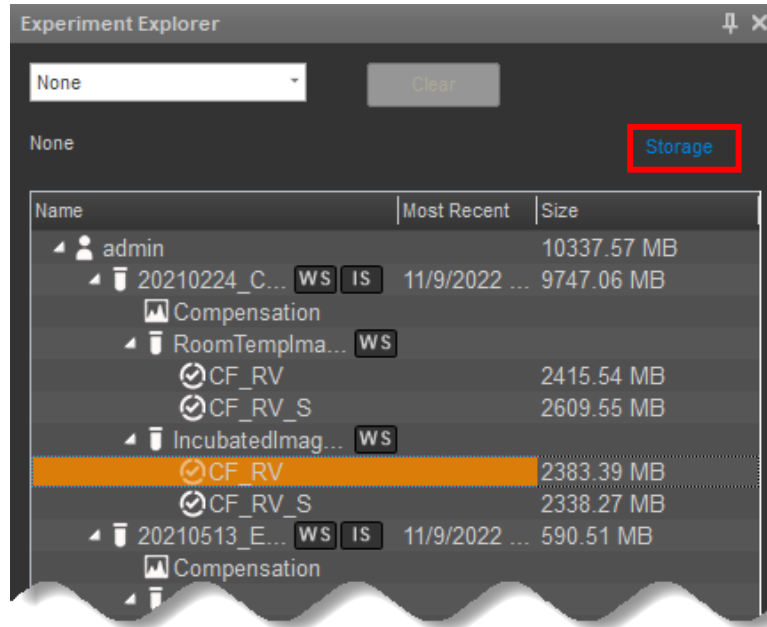


Name	Most Recent	Size
admin		10337.62 MB
20210224_CA... WS IS	11/9/2022 ...	9747.06 MB
Compensation		
RoomTemplmag... WS		
CF_RV		2415.54 MB
CF_RV_S		2609.55 MB
IncubatedImage... WS		
CF_RV		2383.39 MB
CF_RV_S		2338.27 MB
20210513_Ec... WS IS	11/9/2022 ...	590.51 MB
Compensation		
Group		
EcoliCount IS		63.83 MB
EcoliCount(1)		63.82 MB
SepiCount		462.76 MB
Group(1)		
Experiment WS IS CS	2/7/2023 5...	0.03 MB
Plate	2/7/2023 5...	0.03 MB

- By default, the **User** folder is displayed on top of the files list and bears the name of the user currently logged in.
- The files in the folders are arranged in a tree structure, which displays the hierarchy of Plates, Experiments, Groups, and Samples.  
The tree structure varies for Plate and Tube-based Experiments (see “Experiment hierarchy” on page 291).
- Badges next to individual nodes or files in the hierarchy indicate the presence of Workspaces (WS), Instrument Settings (IS), and Compensation Settings (CS).
- The Size column shows the size data for the Experiment and includes the summed size of all Sample data and any Experiment metadata.  
To hide or show the size column, right-click the Experiment Explorer heading row, then deselect or select Size.
- You can use the **folder context menu** (“User and Shared folder context menus” on page 302) to add a new Plate or Tube Experiment to any folder.
- When you create a new Tube or Plate Experiment, the Experiment hierarchy is expanded by default.
- When items are selected in the Experiment Explorer, the entire row is highlighted as specified.
- To sort the Experiments by name, date, or size, click the appropriate column header.

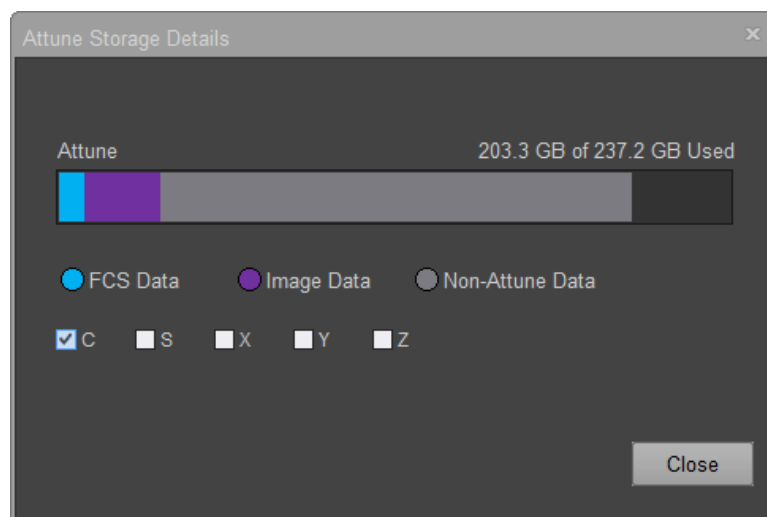
## Storage and Size

To view the data storage details on the instrument, click the **Storage** hyperlink at the top of the **Experiment Explorer**, which opens the **Attune Storage Details** window.

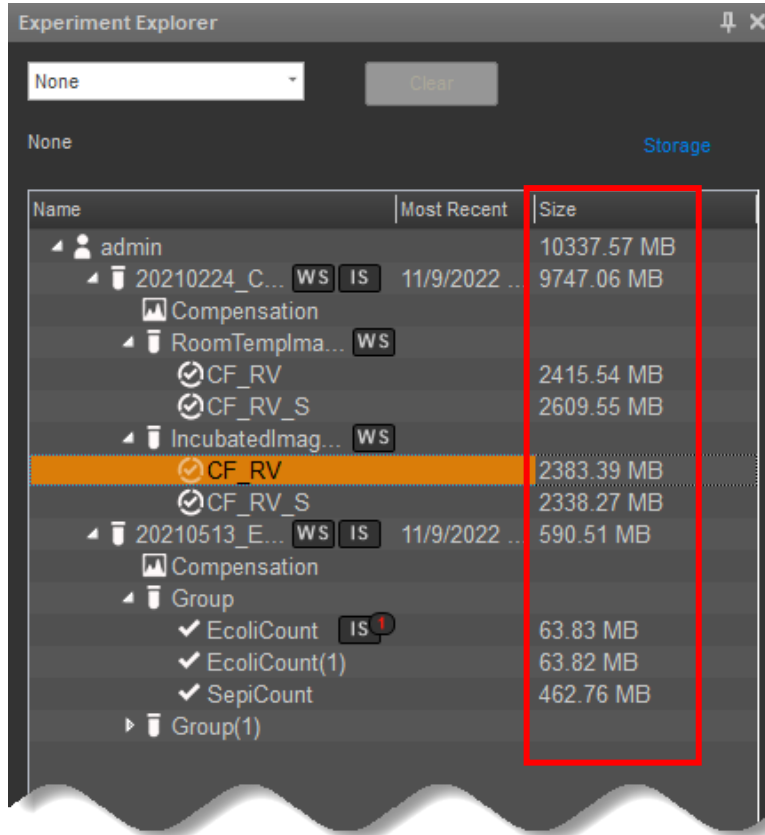


The **Attune Storage Details** displays the total used space and the total combined space broken into:

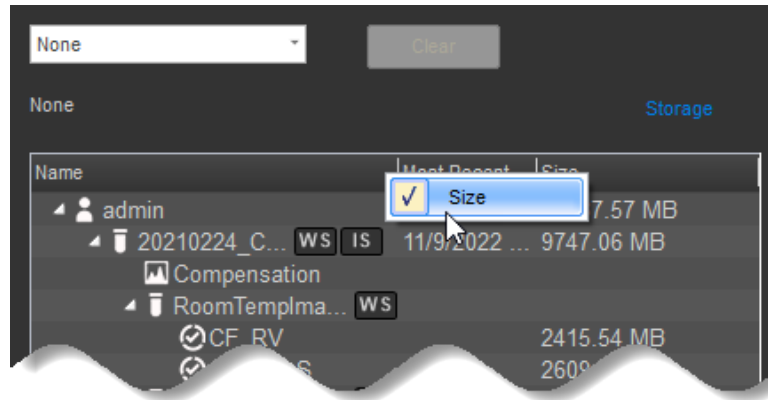
- FCS Data
- Image Data
- Non-Attune Data



The size details for Experiments, Groups, and Samples are shown in the **Size** column in **Experiment Explorer**, which are calculated on sign in and when any changes are made to an experiment (data added or removed).



- The size data are displayed for all samples that contain data and include the size of the FCS file, the image data, and all metadata for the sample.
- The size data are displayed for the Experiment and include the summed size of all Sample data and any Experiment metadata.
- To sort the size data in ascending or descending order, click the **Size column header**.
- To hide or show the size column, right-click the **Experiment Explorer heading row**, then deselect or select **Size**.

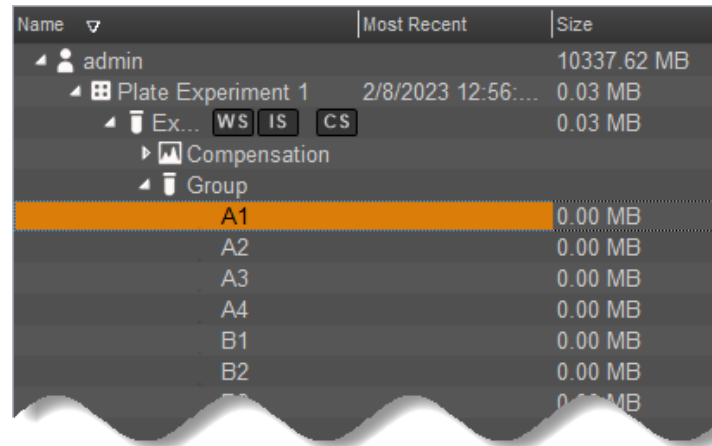




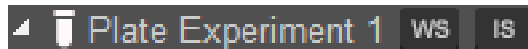
# Experiment hierarchy

## Plate hierarchy

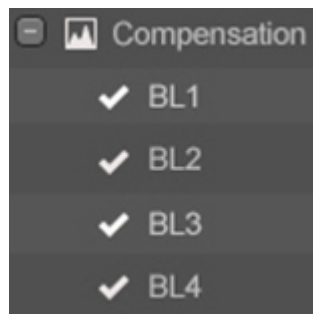
The Experiment Explorer displays the Plate Experiments in a defined hierarchy with a tree structure.



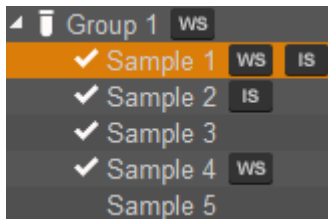
- When you create a new Plate, the software automatically creates a Plate Experiment containing a Compensation node. However, the software does not create any Groups or Samples for new Experiments, unless specified in the *New Experiment dialog* (“New Experiment dialog” on page 606).
- A Plate can contain one Plate Experiment, which has an associated Workspace, as well as Instrument and Compensation settings.  
Badges to the right of the Experiment name indicate the presence of Experiment-level Workspace (WS), Instrument (IS), and Compensation (CS) settings.



- The Experiment-level compensation settings can have multiple associated Compensation samples, which are displayed in laser and detector order.
- When compensation is recorded for a Compensation sample, the check mark icon appears to the left of the Compensation sample name, indicating that the Sample has recorded data.  
If Compensation has not been recorded for a Sample, then no icon appears.



- Each Experiment that has assigned Samples consists of one or more Groups. A Workspace badge to the right of the Group name indicates the presence of Group-level Workspace. Each group can contain up to 400 Samples. Samples can be derived from a well in a Plate or a Tube.



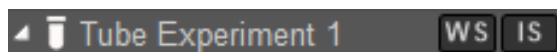
- Each Sample can have a Sample-level Workspace. A Sample can also have Sample-level Instrument settings used instead of the Experiment defaults. Badges to the right of the Sample name indicate the presence of Sample-level Workspace (WS) and Instrument (IS) settings.
- When the Sample is recorded, the check mark icon appears to the left of the Sample name, indicating that the Sample has recorded data. If the Sample has not been recorded, then no icon appears.

## Tube hierarchy

The Experiment Explorer displays the Tube Experiments in a defined hierarchy with a tree structure.



- When you create a new Tube Experiment, the software automatically creates a Compensation node in the hierarchy. However, the software does not create any Groups or Samples for new Experiments, unless specified in the *New Experiment dialog* (“New Experiment dialog” on page 606).
- Each Tube Experiment has an associated Workspace, as well as Instrument and Compensation settings. Badges to the right of the Experiment name indicate the presence of Experiment-level Workspace (WS), Instrument (IS), and Compensation (CS) settings.

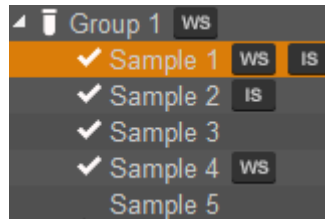


- The Experiment-level compensation settings can have multiple associated Compensation samples, which are displayed in laser and detector order.

- When compensation is recorded for a Compensation sample, the check mark icon appears to the left of the Compensation sample name, indicating that the Sample has recorded data. If compensation has not been recorded for a Sample, then no icon appears.



- Each Experiment that has assigned Samples consists of one or more Groups. A Workspace badge to the right of the Group name indicates the presence of Group-level Workspace. Each group can contain up to 400 Samples. Samples can be derived from a well in a Plate or from a Tube.
- Each Sample can have a Sample-level Workspace. A Sample can also use Sample-level Instrument settings instead of the Experiment defaults. Badges to the right of the Sample name indicate the presence of Sample-level Workspace (WS) and Instrument (IS) settings.







- When the Sample is recorded, the check mark icon appears to the left of the Sample name, indicating that the Sample has recorded data. If the Sample has not been recorded, then no icon appears.

## Experiment Explorer behavior

### Selection indicators

Colors are used to indicate various states of selection in the Experiment Explorer. When items are selected, the entire selected row in the Files view, including the name and date columns, is highlighted.

Selection	Indicator color	Description
Mouse over		When the mouse is moved over an item in the Experiment Explorer, the item is highlighted in gray.
Selected item		When an item is selected in the Experiment Explorer, the item is highlighted in dark blue.
		You can select multiple items simultaneously by holding the <b>Ctrl</b> or the <b>Shift</b> key when clicking. Items which are multiselected are highlighted in light blue, except the last selected item, which is highlighted in dark blue.
Active item		When an item is loaded into the current Workspace, the Sample name in Experiment Explorer is highlighted in orange.

### Selection

- Double-click the **Plate node** or **Plate** or **Tube Experiment node** to set the focus to the **Heat Map view** (Chapter 6, “Heat Map View”) for that Experiment.  
If another Experiment was already open, this action opens the first Sample in the selected Experiment.  
If no Sample exists in the selected Experiment (if Plate), then double-clicking the node opens the **Experiment-level Workspace** and sets the **Experiment node** as active. The **Workspace dropdown** is disabled.
- Double-click a **Group node** to activate the first Sample in the Group.  
If another Experiment is already open, this action opens the first Sample in the selected Group.  
If no Sample exists in the selected Group, then double-clicking the node opens the **Experiment-level Workspace** and sets the **Experiment node** as active. The **Workspace dropdown** is disabled.
- Double-click the **Compensation node** or any **Compensation sample** to set the focus to the **Compensation Workspace** for the selected Sample. The **Compensation tab** becomes available in the Ribbon bar (“Experiment tab controls” on page 95).
- If you double-click the main **Compensation node** and Compensation controls exist, the first Compensation control Sample is selected and the focus is set to the **Workspace view** (Chapter 5, “Workspace view”).  
If no Compensation controls exist, double-clicking the Compensation node opens the **Compensation Setup dialog** described on “Compensation setup dialog” on page 581.  
Double-clicking the **Compensation node** that shows the “read-only” icon (🔒) opens the **Compensation Matrix** (“Matrix dialog” on page 599) for that Experiment.

- Double-clicking a **Sample** opens the Sample using the currently selected Workspace and loads the associated data, if present.  
If a Sample does not have the selected Workspace, it reverts to the next Workspace in the hierarchy (Sample→Group, or Group→Experiment).  
If a new Experiment is loaded, the Workspace is set to the Experiment-level Workspace.
- During acquisition, the active Sample cannot be changed
- During acquisition, double-clicking the active **Sample's Plate node**, or the **Plate** or **Tube Experiment node** sets the focus to the **Heat Map view** for the active Experiment.
- The double-clicking any other node or object does not have an effect.

## Drag and drop

### Workspace (WS)

A Workspace (WS) badge can be dragged from an Experiment, Group, or Sample, and dropped to another Experiment, Group, or Sample.

- To apply the Workspace from one location (Experiment, Group, or Sample) to another location, click the **WS badge**, then drag the badge to the destination location while continuing to hold on the badge. Release the WS badge at the new location to apply the Workspace to the location.
- Dragging and dropping a WS badge in the Experiment Explorer creates a copy of the Workspace at the dropped location. The copy of the Workspace has contents identical to the original, but it is not linked to it.
- Experiment-level WS badge can be dragged to another Experiment, Group, or Sample, which copies the Workspace to the Experiment, Group, or Sample.
- Group-level WS badge can be dragged to another Experiment, Group, or Sample, which copies the Workspace to that respective level.
- A Sample-level WS badge can be dragged to another Experiment, Group, or Sample, which copies the Workspace to that respective level.

### Instrument Settings (IS)

An Instrument Settings (IS) badge can be dragged from an Experiment or Sample to another Experiment, Group, or Sample.

- To apply the Instruments Settings from one location (Experiment, Group, or Sample) to another location, click the **IS badge**, then drag the badge to the destination location while continuing to hold on the badge. Release the IS badge at the new location to apply the instrument settings to the new location.
- Experiment-level IS badge can be dragged to another Experiment or empty Samples, which updates the IS for the Experiment or the Samples.  
If the Experiment contains acquired Samples that have Experiment-level IS, the Samples are updated to show the Sample IS indicator.
- Experiment-level IS badge can also be dragged to a Group folder to apply the Experiment IS to all unrecorded Samples in that Group.  
If this is done within the same Experiment, any Sample-level IS that previously existed in that Group is cleared.
- Experiment-level IS badge cannot be dragged to an Experiment that has existing compensation.

- If the Experiment contains acquired Samples that already have Sample-level IS, dragging the IS badge has no effect on these Samples.
- Sample-level IS badge can be dragged to other Experiments or empty Samples.
- Sample-level IS badge can be dragged to a Group folder to apply the IS to all unrecorded Samples in that Group.
- Dragging a Sample-level IS badge to an Experiment updates the Experiment IS. When this is done, the Sample-level IS badge is removed from the Sample, which is set to use the Experiment-level IS. All other Sample-level IS badges remain regardless of their similarity to the current Experiment-level IS.
- Sample-level IS badge cannot be dragged to a Sample with a recorded FCS file.
- The IS badge cannot be dropped on an Experiment, Group, or Sample, where the instrument configuration does not match. In this case, a warning dialog is displayed with the message: “The instrument settings are not supported by the current instrument configuration and cannot be imported”.

## Compensation Settings (CS)

A Compensation Settings (CS) badge can be dragged from one Experiment to another Experiment.

- To apply the Compensation Settings from one Experiment to another, click the **CS badge**, then drag the badge to the destination Experiment in the Experiment Explorer while continuing to hold on the CS badge. Release the CS badge at the new location to apply the Compensation and Instrument settings from the original Experiment to the new Experiment.
- The CS badge is only shown next to an Experiment, if the Experiment has compensation defined.
- The CS badge can be dragged from one Experiment to another, which copies the Compensation Settings and the corresponding Instrument Settings to that Experiment.
- When Compensation Settings are applied to another Experiment via drag-and-drop, they are applied as read-only compensation based on an XML file. This is indicated in the software by a padlock icon next to the Compensation Control.
- The original FCS files recorded for the Compensation Controls are not viewable from the Experiment in which the settings were copied; only the compensation values are visible in the Compensation Matrix (“Matrix dialog” on page 599)
- The Compensation Settings can only be applied to another Experiment, if the instrument models match and the target Experiment does not already have compensation defined or recorded.
- The rules for drag-and-drop of the Compensation Settings badge follow the same guidelines described for dragging an IS badge to an Experiment-level IS.

## Instrument Settings (IS) badge identification

The **Instrument Settings (IS) badge** is used to identify Samples that have Sample-level Instrument Settings. To provide extra information whether the Samples in an Experiment have different Instrument Settings, a number is placed with the IS badge to show which Samples have the same Instrument Settings.

- When samples show an IS badge, they are compared with each other to determine whether the Instrument Settings XML is different.
- If the Sample-level IS differs from the Experiment-level IS, the Sample IS displays an extra badge with a colored number to indicate that the Instrument Settings are different (for example, IS1, IS2, etc.).



- If the Instrument Settings are the same as the experiment level IS, the IS badge is displayed without the extra colored number badge.
- Each Sample is compared to the Experiment-level Instrument Settings and to the other Sample Instrument Settings (this also applies when pasting IS settings via the Heat Map or when dragging and dropping the IS badges in the Experiment Explorer).
- Each Sample with new or unique Instrument Settings is assigned an IS badge that includes a new number (base 1), which is incremented sequentially. All Instrument Settings that are shared show the same colored number badge.

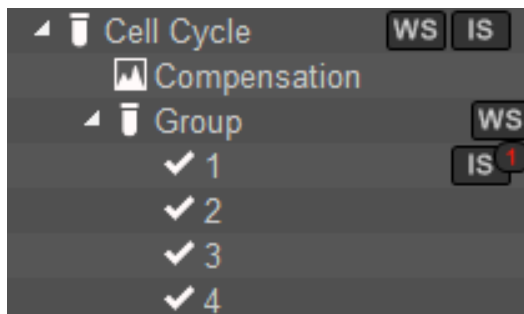


Figure 49 Sample 1 has a unique, Sample-level IS that differs from the Experiment-level IS.

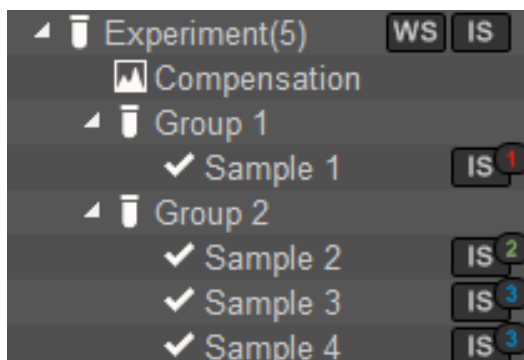


Figure 50 Samples 1, 2, and 3 have unique ISs

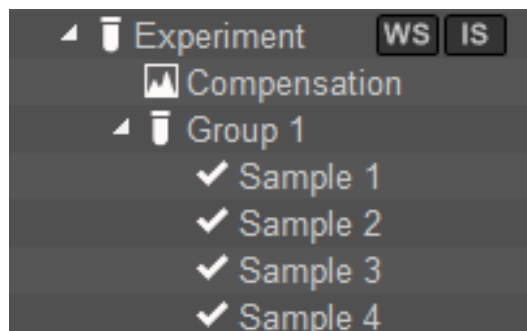


Figure 51 Samples 1, 2, 3, 4 have the same IS

- If the Sample-level IS badge with a number is dragged to Experiment-level, the Sample-level IS number is removed so that only the IS badge is displayed (that Sample IS now equals the Experiment-level IS), and the remaining badges are renumbered.

## Drag-and-drop from Experiment Explorer to Overlays

You can select Samples to drag-and-drop on to plots in the Overlay gallery view (Chapter 9, “Overlays”).

- When Samples are dragged from the Experiment Explorer to a plot in the Overlay gallery view, new plots are added to the gallery corresponding to the selected Samples and plot.
- The plot into which the dragged Samples are dropped is used to set the gallery plot customization for the newly added plots.

## Drag-and-drop FCS files

You can import a Sample by dragging an FCS file onto an empty Sample in the Experiment Explorer.

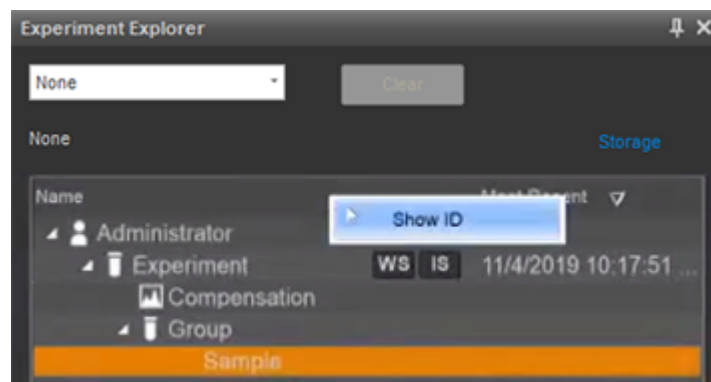
- If the drag-and-drop operation is valid, the drop icon (a + sign on top of the arrow) is displayed.
- If the Sample already has data, Acquisition is in progress, Automation mode is enabled, or the instrument is paused, the drop option is disabled and the drop icon is displayed (the no entry icon is displayed).
- When the FCS file is dropped onto the empty Experiment Explorer item, the FCS file is imported as though it was imported using the Import FCS File command in the Sample context menu (“Import FCS file” on page 326). The Sample name in the Experiment Explorer is updated to the name of the FCS file after the import.
- Only a valid FCS file can be imported.



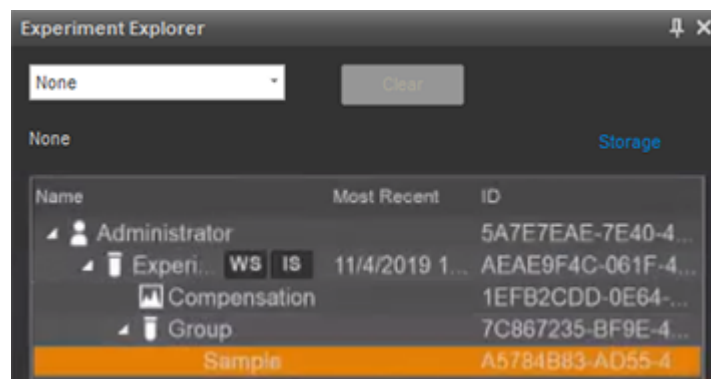
## Display ID

When the Attune™ Cytometric Software is in the SAE mode, you can show the ID of application objects in the Experiment Explorer.

- Application objects are the domain objects that are of interest for auditing and e-signing when the Attune™ Cytometric Software is in the SAE mode (see Appendix D, “SAE Administrator Console”).
- Examples of application types in the Attune™ Cytometric Software are Plate Experiment, Tube Experiment, Template, Filter Configuration, and Filter Definition.
- The GUID is the unique ID assigned each an application object. It is an immutable property and cannot be changed.
- To show the GUID of application objects in Experiment Explorer, right-click the column header in Experiment Explorer, then select **Show ID**.



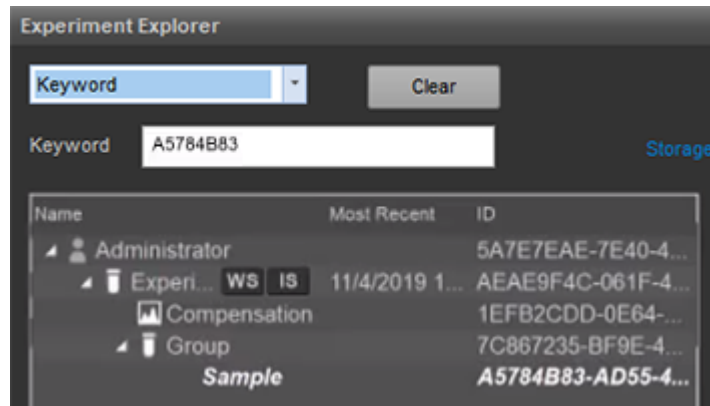
The GUID for the application objects are displayed in the ID column of Experiment Explorer.



- When the ID column is shown on the Experiment Explorer, you can use the GUID to search for specific application objects.

To do this, select **Keyword** in the Search dropdown, enter the GUID of the application object you want to search, then click **Enter**.

If found, the application object you have searched appears bolded and italicized in the Experiment Explorer.



# Experiment Explorer context menus

## Overview

Experiment Explorer has context menus associated with different levels of the hierarchy in the tree structure. You can access each context menu by right-clicking the appropriate object in the tree structure.

- The Experiment Explorer context menus include the following:
  - *User context menu* (“User and Shared folder context menus” on page 302)
  - *Plate context menu* (“Plate context menu” on page 304)
  - *Experiment context menu* (“Experiment context menu” on page 310)
  - *Group context menu* (“Group context menu” on page 318)
  - *Sample context menu* (“Sample context menu” on page 324)
  - *Compensation node context menu* (“Compensation node context menu” on page 331)
  - *Compensation control context menu* (“Compensation control context menu” on page 335)
- Unless indicated otherwise, all Experiment Explorer context menu items are always visible and enabled.
- When a combination of different objects is selected in the Experiment Explorer (e.g. Experiment, Group, and Samples), the right-click context menus are not available.

## Validation of name fields

Context menu options that allow you to create or edit names of items follow the validation rules outlined below:

- **Duplicate** context menu option duplicates the selected object and displays the name in a textbox in the Edit mode.
- **Rename** context menu option displays the current name of the selected object in a textbox in the Edit mode.
- You can rename a node by pressing the **F2** key to enter the Edit mode.
- Click anywhere outside the edit textbox or press **Enter** to validate the entered name and close the Edit mode.
- If a valid name is present in the textbox, the name is updated.
- Leading and trailing spaces are removed on validation. Consecutive spaces are converted to single spaces on validation.
- If you attempt to enter invalid characters into the edit textbox, a warning dialog indicates the error condition and the invalid characters do not appear in the textbox.
- If an invalid name is present when focus is lost, then the name reverts to the previous valid name.
- If the entered name already exists, when focus is lost, a warning dialog indicates that the entered name already exists and the textbox remains in Edit mode.

## User and Shared folder context menus

Right-click the **User folder heading** to show the **User folder** or the **Shared folder context menu**, which enable you to create New Experiments from scratch or using a Template, or to import predefined Experiments.

During acquisition, only the **Collapse All** option is enabled.

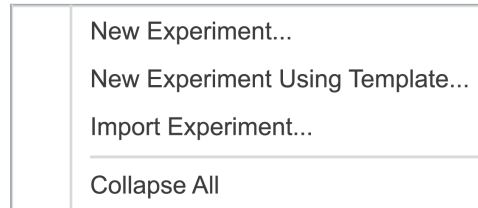


Figure 52 User folder context menu

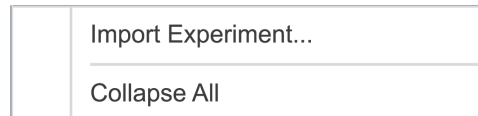


Figure 53 Shared folder context menu

### New Experiment

- **New Experiment** opens the **New Experiment dialog** (“New Experiment dialog” on page 606), which enables you to create a new Plate or Tube Experiment or to import FCS files into a new Analysis Experiment.
- The **New Experiment** option is not available on any Shared folder context menus. It is disabled if a Plate is paused or if the application is in the Automation mode.

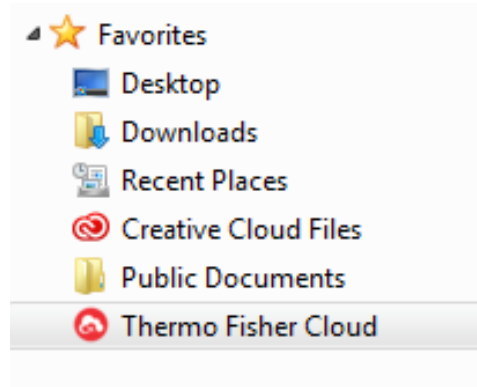
### New Experiment Using Template

- **New Experiment Using Template** opens the **New Experiment from Template dialog** (“New experiment from template dialog” on page 613), which enables you to select a Plate- or Tube-based Experiment as a template for a new Experiment.
- The **New Experiment Using Template** option is disabled if a Plate is paused or if the application is in the Automation mode.

### Import Experiment

- **Import Experiment** opens the **File Open (Import) dialog** (“File Open (Import) dialog” on page 721), which enables you to import preexisting Plate- or Tube-based Experiments.
- The **Import Experiment** option is not available during acquisition.

- If the device (instrument or software) has been registered to a Connect account, the File Open dialogue includes an icon for the Connect listed under Favorites. Clicking on this icon opens a virtual folder listing all files and folders included in the Connect account. You can import the Experiment from the Connect account by selecting the desired location in the Connect account during Experiment import.



- Connect account contents are only visible when viewed through the **File Open (Import)** and **File Save (Export)** dialogues from the Attune™ Cytometric Software.

## Collapse All

- **Collapse All** collapses all expanded nodes down to the Plate or Tube Experiment-levels for the current Folder.

Name	Most Recent	Size
admin		10337.65 MB
▶ Tube Experim...	WS IS 2/8/2023 ...	0.03 MB
▶ Plate Experiment 1	2/8/2023 ...	0.03 MB
▶ Experiment	WS IS CS 2/7/2023 ...	0.03 MB
▶ 20210513_Ec...	WS IS 11/9/2022 ...	590.51 MB
▶ 20210224_CA...	WS IS 11/9/2022 ...	9747.06 MB

The **Collapse All** option is always available.

## Plate context menu

Right-click a **Plate icon** or **Plate name** in **Experiment Explorer** to show the **Plate context menu**.

During acquisition, all options in **Plate context menus** are disabled.

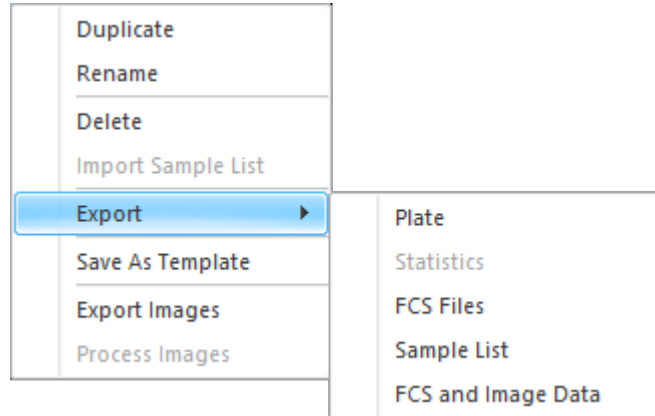


Figure 54 Plate context menu in User folder

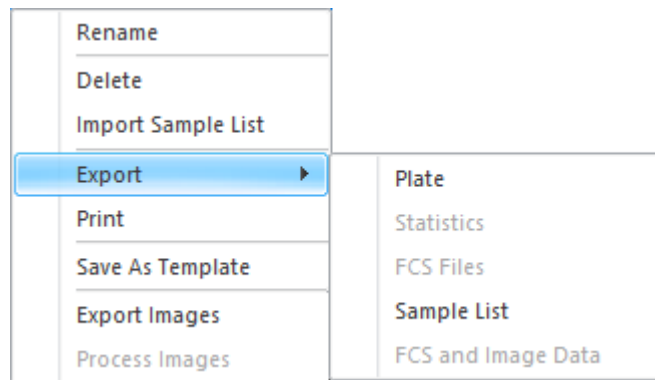


Figure 55 Plate context menu in Shared folder

## Duplicate

- **Duplicate** enables you to duplicate the selected plates and their full hierarchy (i.e., Experiment, Compensation, Groups, and Samples).
- The **Duplicate** option is enabled when single or multiple plates are selected. This option is disabled if the Plate Experiment Instrument settings do not match the system configuration
- When a Plate Experiment is duplicated, all Workspaces (Experiment-, Group-, and Sample-level), Run Protocols, and annotations are preserved and all Experiment and Sample-level keywords are duplicated. Only the Experiment-level Instrument Settings are preserved. Sample-level Instrument settings and FCS files are not duplicated.
- If Compensation settings are available, the newly created Plate contains Experiments with empty Compensation wells or Tubes as defined in the source Experiment.

- The newly created Plate must have a unique name. When a Plate is duplicated, it is automatically assigned a name, which consists of the same name as the current Plate plus an incrementing numerical suffix. The numerical suffix starts at 1 and increments until a unique identifier is achieved.
- The Plate name appears in an edit textbox ready for editing. The newly entered name is validated as described on “Validation of name fields” on page 301.

## Rename

- **Rename** opens an edit textbox containing the name of the current Plate in the edit mode.
- The **Rename** option is enabled when a single plate is selected.
- Clicking anywhere outside the edit textbox validates the Plate name as described on “Validation of name fields” on page 301.
- You can also rename a selected Plate by pressing the **F2** key on the keyboard to enter the editing mode.

## Delete

- **Delete** opens the **Deleted Items dialog** (“Deleted items dialog” on page 339), which enables you to delete the selected Plates.
- The **Delete** option is enabled when single or multiple Plates are selected.

## Import Sample List

**Import Sample List** opens the **File Open (Import) dialog** (“File Open (Import) dialog” on page 721) with the extension set as CSV by default (\*.csv), which enables you to import a Sample List from an existing CSV file. Other file formats are not supported.

- The **Import Sample List** option is disabled if the Experiment contains any recorded data, multiple nodes of the same level are selected in Experiment Explorer, acquisition is in progress, or the Automation mode is enabled.

When the Sample List is imported, the software looks for the following columns in the CSV file:

LOCATION	EXPERIMENT
SAMPLE	EXP_NOTES
SAMPLE_NOTES	PLATE
GROUP	PLATE_ID
GROUP_NOTES	PLATE_NOTES

LOCATION column is required for import and must contain data in the format A1, A2, A3,... etc. (for a Plate Experiment) or T1, T2, T3,... etc. (for a Tube Experiment). All other columns are optional.

PLATE, PLATE\_ID, and PLATE\_NOTES columns are only created when importing into a Plate Experiment. These columns are ignored when importing the Sample List into a Tube Experiment.

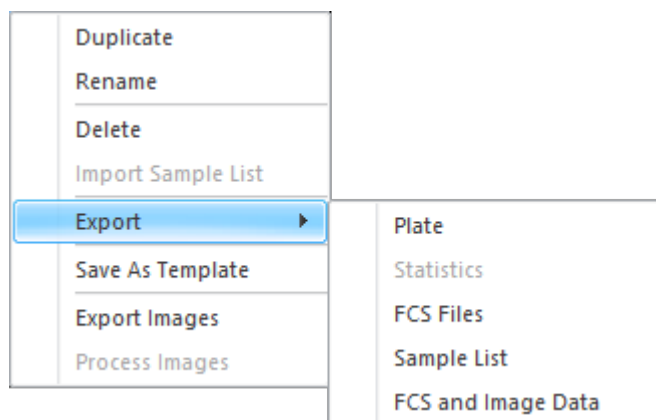
- When importing the CSV sample list, the **Map Sample List dialog** (“Map Sample List Data dialog” on page 764) is displayed, which enables you to create and map Samples based on Location (i.e., a Sample mapped to A1 in the Sample List will be mapped to A1 on the Heat Map).

Extra keywords are mapped based on the keywords defined for the Experiment. If the Experiment does not have the custom keywords included in the Sample List CSV file, the custom keywords in the file are not added to the Sample List in the Experiment.

- If Samples already exist in the Experiment or Plate (i.e., Samples are assigned to a location), a prompt to update the Sample information is displayed (“Update sample information dialog” on page 766).
- After the existing Samples are updated, the import process continues to process the addition of new Samples, as necessary.
- The Sample names created in each Group during the import process must be unique, and Sample, Group, Experiment, and Plate names must be less than or equal to 50 characters long.
- Names cannot contain the illegal characters `/:*? "<>|&` and end with a period.  
Names also cannot be any of the following words: CON, PRN, AUX, CLOCK\$, NUL, COM1, COM2, COM3, COM4, COM5, COM6, COM7, COM8, COM9, LPT1, LPT2, LPT3, LPT4, LPT5, LPT6, LPT7, LPT8, and LPT9.

## Export

**Export** option displays a submenu containing **Plate**, **Statistics**, **FCS Files**, **Sample List**, and **FCS and Image Data** as options. The **Export** option is enabled when single or multiple Plates are selected.



- **Export ▶ Plate** opens the **File Save (Export) dialog** (“File Save (Export) dialog” on page 715), which enables you to save the selected Plates in the desired location.
- **Export ▶ Statistics** opens the **Export Statistics dialog** (“Export Statistics dialog” on page 737), which enables you to export the statistics of selected samples and selected views as a single file or as individual files.
  - Statistics for all Samples in the selected Experiment are exported except Samples with no data.
  - Statistics can be exported as a single file (statistics for all Samples are combined into a single CSV file) or as individual files (statistics for all Samples are exported into separate CSV files).
  - The **Export ▶ Statistics** option is always enabled, except during acquisition.



- **Export ▶ FCS Files** opens the **File Save (Export) dialog**, which enables you to save the selected FCS files using the required FCS version in the desired location.
  - All FCS files in the selected Experiments are exported. If a Sample in the selected Experiment has no FCS file, then no exported FCS file is created for that Sample.
  - If the Experiment compensation settings are different from the Compensation settings contained in the FCS file, the target and label names have been changed, or Experiment or custom keywords have been changed, a prompt to update the keywords is displayed.
  - The **Export ▶ FCS Files** option is disabled during data acquisition (Run or Record) or if the selected Experiment does not contain FCS data files.
- **Export ▶ Sample List** opens the **File Save (Export) dialog**, which enables you to save the Sample List in the desired location.
  - The **Export ▶ Sample List** option is disabled if acquisition is in progress or the Automation mode is enabled. In both cases, this option is enabled when a plate is paused. If multiple nodes of the same level are selected, this option is also disabled.
  - When the Sample List is exported, a CSV file is created that contains a list of all Samples in the Experiment (Compensation controls are not included in the Sample List) with the following columns:

LOCATION	EXPERIMENT
SAMPLE	EXP_NOTES
SAMPLE_NOTES	PLATE
GROUP	PLATE_ID
GROUP_NOTES	PLATE_NOTES

- PLATE, PLATE\_ID, and PLATE\_NOTES columns are only created when exporting from a Plate or Plate Experiment. These columns are not created when exporting the Sample List from a Tube Experiment.
  - All user created keywords are also included in the export set, and the user defined keyword names are used as the column headers for the user defined keywords.
  - If a field is blank, it is left empty in the resulting CSV file.
  - For Tube samples, the location information is based on Tube position in the Heat Map and is designated as T1, T2, etc.
- **Export ▶ FCS and Image Data** opens the **File Save (Export) dialog**, which enables you to save the FCS files and imaging data together in the desired location.
    - By default, images, extended parameters (image processing data), and image mask are exported into a single zip file based on the ACS file format.
    - Exported ACS files can be imported into a sample, which imports the FCS data and all imaging data.
    - Exported ACS files can be opened in third party software that supports the ACS standard (such as FlowJo™ Software and FCS Express™ Software).

---

**IMPORTANT!** The **Export ▶ FCS and Image Data** option is available only for experiments performed with an Attune™ CytPix™ Flow Cytometer, and it is active only if the experiment has FCS and image data.

---

## Save as Template

**Save As Template** opens the **Save As Template dialog** (“Save As Template dialog” on page 747), which enables you to name the current Plate and save it as a template in database for future reuse.

## Export Images

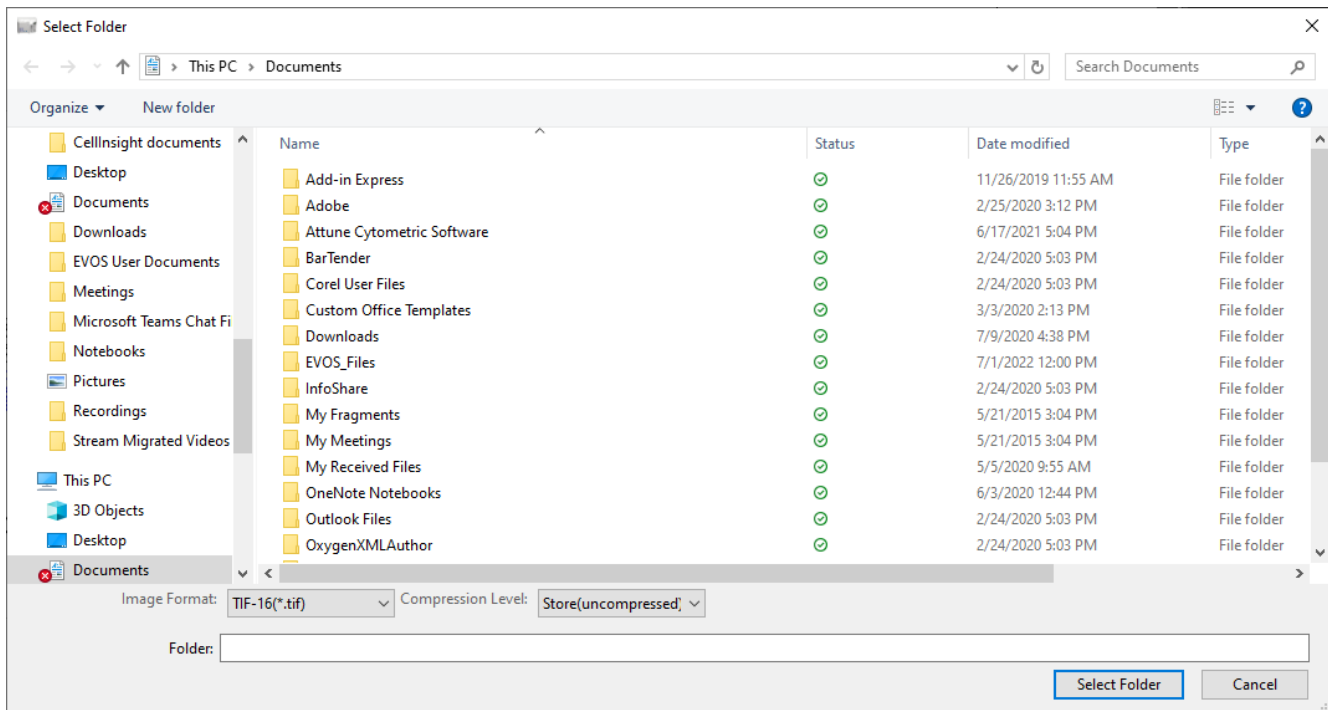
**Export Images** opens the **Select Folder (Folder browser) dialog** (“Folder Browser dialog” on page 723), which enables you to save the images from the selected Plate in the desired location.

**Note:** The **Export Images** option is available only for experiments performed with an Attune™ CytPix™ Flow Cytometer, and it is active only if the experiment has image data.

- When images are exported, all images for the selected sample or samples are exported into a single zip file.
- By default, images are exported as **TIF-16** files, which is the native file format collected from the Attune™ CytPix™ Flow Cytometer.
- You can select to save the images as **TIF-16, TIF-8, PNG, BMP, GIF, EMF, or JPG** files using the **Image Format** dropdown in the **Select Folder (Folder browser) dialog**.

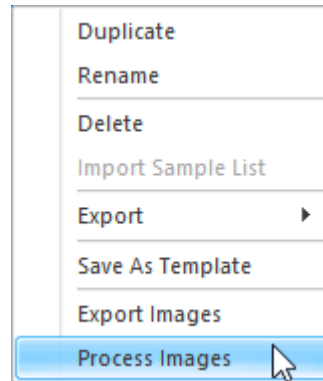
**IMPORTANT!** When saving in any format other than **TIF-16**, images are converted to 8-bit images, which results in loss of image pixel data.

- Use the **Compression Level** dropdown to select to store the images uncompressed (**Store (uncompressed)**) or save them **Compressed** to save disk storage space.



## Process Images

**Process Images** opens the **Process Images** dialog (see Chapter 23, “Process Images dialog”), which enables you to process captured images from selected **Samples** and **Populations** using system-provided **Image Processing Models**.



- The **Process Images** option is only visible for Attune™ CytPix™ Experiments.
- You can only process images for the Experiment that is active in the Workspace. The **Process Images** option is greyed out in the **Plate context menus** of Experiments that are not active.

---

**Note:** For more information about the Attune™ Cytometric Software image processing workflow, see the *Attune™ Cytometric Software Image Processing Workflow* (Pub. No. [MAN0028531](#)), which is available for download at [thermofisher.com](#).

---

## Experiment context menu

- Right-click a **Tube Experiment node** in **Experiment Explorer** to open the **Tube Experiment context menu**.
- Right-click a **Plate Experiment node** in **Experiment Explorer** to open the **Plate Experiment context menu**.
- During acquisition, all options in **Experiment context menus** are disabled.

**Note:** The **Export Images** and **Process Images** options are available only for experiments performed with the Attune™ CytPix™ Flow Cytometer. These options are not visible in Attune™ NxT experiments.

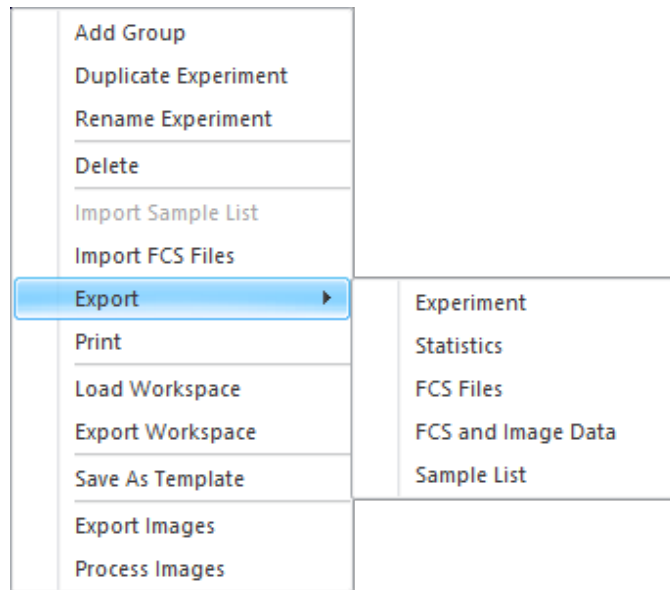


Figure 56 Tube Experiment context menu

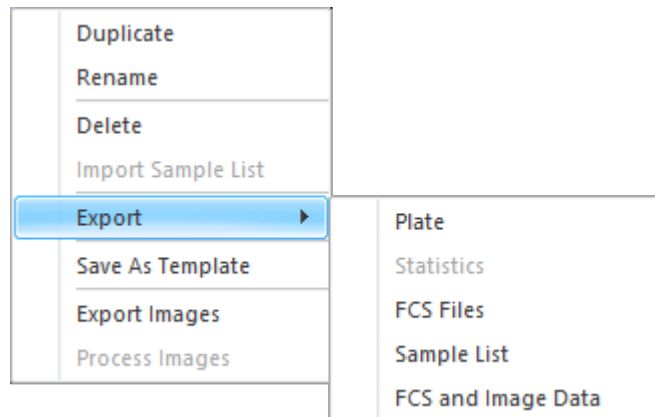


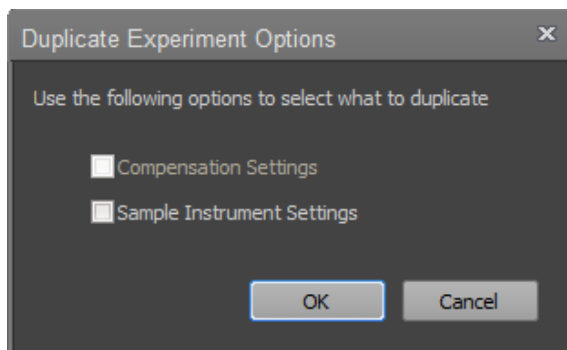
Figure 57 Plate Experiment context menu

## Add Group

- **Add Group** opens the **New Group dialog** (“New experiment from template dialog” on page 613), which enables you to add a Group to the end of the currently selected Experiment.
- You can create up to 100 Groups for each Experiment.
- **Add Group** option is only enabled when a single Tube Experiment is selected.

## Duplicate Experiment

- **Duplicate Experiment** duplicates the selected Experiments with their associated Groups and Samples.
- The **Duplicate Experiment** option is visible only for Tube Experiments, and it is enabled only when single or multiple Experiments are selected.  
This option is disabled if the Experiment Instrument settings do not match the system configuration
- When an Experiment is duplicated, all Workspaces (Experiment-, Group-, and Sample-level), Run Protocols, and annotations are preserved.  
Experiment-level Instrument Settings (IS) are preserved. If Sample-level IS exists in the Experiment, a duplicate Experiment options menu provides the option to duplicate Sample-level IS and Compensation Settings.



FCS files are not duplicated.

All Experiment- and Sample-level keywords are duplicated.

- If Compensation Settings have not been duplicated, the newly created Experiment contains a Compensation node with empty Samples as defined in the source Experiment.  
If the Compensation Settings have been duplicated, the newly created Experiment contains a Compensation node with locked Compensation controls matching the compensation values in the source Experiment.
- The newly created Experiment must have a unique name. After an Experiment is duplicated, it is assigned a name that consists of the same name as the current Experiment plus an incrementing numerical suffix. The numerical suffix starts at 1 and increases in increments until a unique identifier is achieved.

## Rename Experiment

- **Rename** opens an edit textbox containing the name of the current Experiment in the edit mode.
- The **Rename** option is enabled when a single Experiment is selected.
- Clicking anywhere outside the edit textbox validates the Experiment name as described on “Validation of name fields” on page 301.
- You can also rename an Experiment by first selecting it, then pressing the **F2** key to enter the editing mode.

## Delete

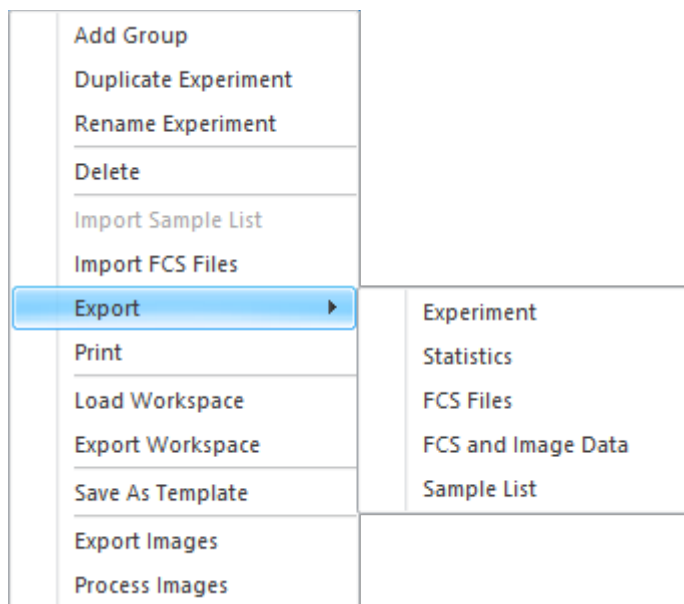
- **Delete** opens the **Deleted Items dialog** (“Deleted items dialog” on page 339), which enables you to delete the selected Experiments.
- The **Delete** option is enabled when single or multiple Experiments are selected.

## Import Sample List

- **Import Sample List** opens the **File Open (Import) dialog** (“File Open (Import) dialog” on page 721) with the extension set as CSV by default (\*.csv), which enables you to import a Sample List from an existing CSV file as described for “Import Sample List” in “Plate context menu” (“Import Sample List” on page 305). File formats other than CSV are not accepted.

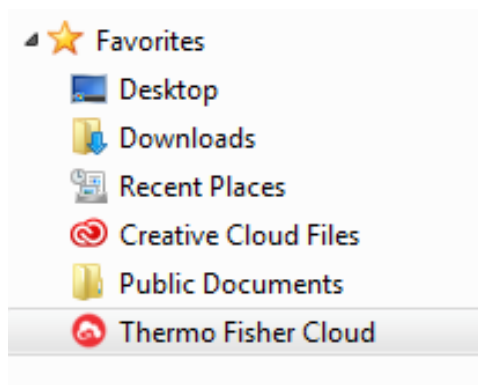
## Export

**Export** displays a submenu containing **Experiment**, **Statistics**, **FCS Files**, **FCS and Image Data**, and **Sample List** as options.

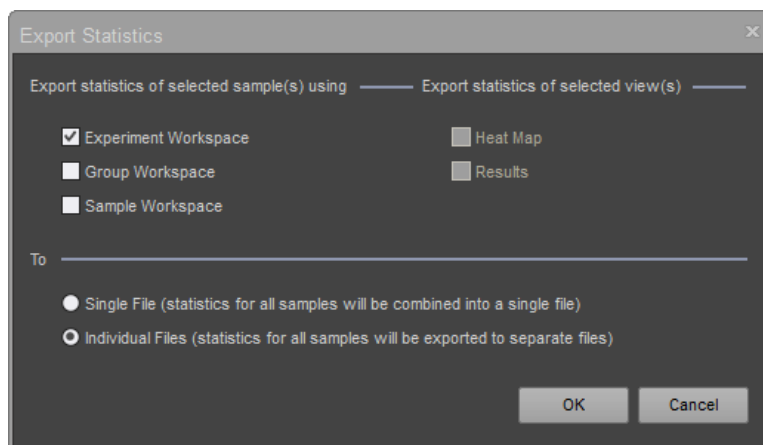


The **Export** option is enabled when single or multiple Plates are selected.

- **Export ▶ Experiment** opens the **File Save (Export) dialog** (“File Save (Export) dialog” on page 715), which enables you to save the selected Experiments in the desired location.
  - The **Export ▶ Experiment** option is visible and enabled only for Tube Experiments.
  - If the device (instrument or software) has been registered to a Connect account, the **File Open dialogue** includes an icon for the Connect listed under **Favorites**. Clicking this icon opens a virtual folder listing all files and folders included in the Connect account. You can save the Experiment to the Connect account by selecting the desired location in the Connect account during Experiment export.



- Connect account contents are only visible when viewed through the **File Open (Import)** and **File Save (Export) dialogues** from the Attune™ Cytometric Software.
- **Export ▶ Statistics** opens the **Export Statistics dialog** (“Export Statistics dialog” on page 737), which enables you to export the statistics of selected samples and selected views as a single file or as individual files.



- Statistics for all Samples in the selected Experiment are exported except Samples with no data.
- Statistics can be exported as a single file (statistics for all Samples are combined into a single CSV file) or as individual files (statistics for all Samples are exported into separate CSV files).
- The **Export ▶ Statistics** option is always enabled, except during acquisition.

- **Export ▶ FCS Files** opens the **File Save (Export) dialog**, which enables you to save the selected FCS files using the required FCS version in the desired location.
  - All FCS files in the selected Experiments are exported. If a Sample in the selected Experiment has no FCS file, then no exported FCS file is created for that Sample.
  - If the Experiment compensation settings are different from the Compensation settings contained in the FCS file, the target and label names have been changed, or Experiment or custom keywords have been changed, a prompt to update the keywords is displayed.
  - The **Export ▶ FCS Files** option is disabled during data acquisition (Run or Record) or if the selected Experiment does not contain FCS data files.
- **Export ▶ FCS and Image Data** opens the **File Save (Export) dialog**, which enables you to save the FCS files and imaging data together in the desired location.
  - By default, images, extended parameters (image processing data), and image mask are exported into a single zip file based on the ACS file format.
  - Exported ACS files can be imported into a sample, which imports the FCS data and all imaging data.
  - Exported ACS files can be opened in third party software that supports the ACS standard (such as FlowJo™ Software and FCS Express™ Software).

---

**IMPORTANT!** The **Export ▶ FCS and Image Data** option is available only for experiments performed with an Attune™ CytPix™ Flow Cytometer, and it is active only if the experiment has FCS and image data.

---

- **Export ▶ Sample List** opens the **File Save (Export) dialog**, which enables you to save the Sample List in the desired location.
  - The **Export ▶ Sample List** option is disabled if acquisition is in progress or the Automation mode is enabled. In both cases, this option is enabled when a plate is paused. If multiple nodes of the same level are selected, this option is also disabled.
  - When the Sample List is exported, a CSV file is created that contains a list of all Samples in the Experiment (Compensation controls are not included in the Sample List) with the following columns:

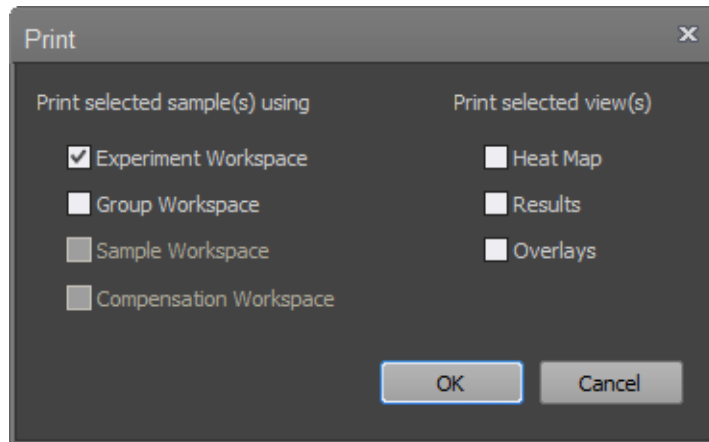
LOCATION	EXPERIMENT
SAMPLE	EXP_NOTES
SAMPLE_NOTES	PLATE
GROUP	PLATE_ID
GROUP_NOTES	PLATE_NOTES

- PLATE, PLATE\_ID, and PLATE\_NOTES columns are only created when exporting from a Plate or Plate Experiment. These columns are not created when exporting the Sample List from a Tube Experiment.
- All user created keywords are also included in the export set, and the user defined keyword names are used as the column headers for the user defined keywords.
- If a field is blank, it is left empty in the resulting CSV file.
- For Tube samples, the location information is based on Tube position in the Heat Map and is designated as T1, T2, etc.



## Print

- **Print** opens the **Batch Print dialog** (“Batch Print dialog” on page 734), which enables you to select a Workspace (Experiment, Group, or Sample) or a View (Heat Map, Results, or Overlays) to print.



- The **Print** option is only enabled when a single Experiment is selected. It is disabled during acquisition.

## Load Workspace

- **Load Workspace** opens the **File Open (Import)** dialog (“File Open (Import) dialog” on page 721), which enables you to select a **Workspace (.aws file)** to import.
- The selected Workspace replaces the existing Experiment-level Workspace.
- The **Load Workspace** option is only enabled when a single Experiment is selected.

## Export Workspace

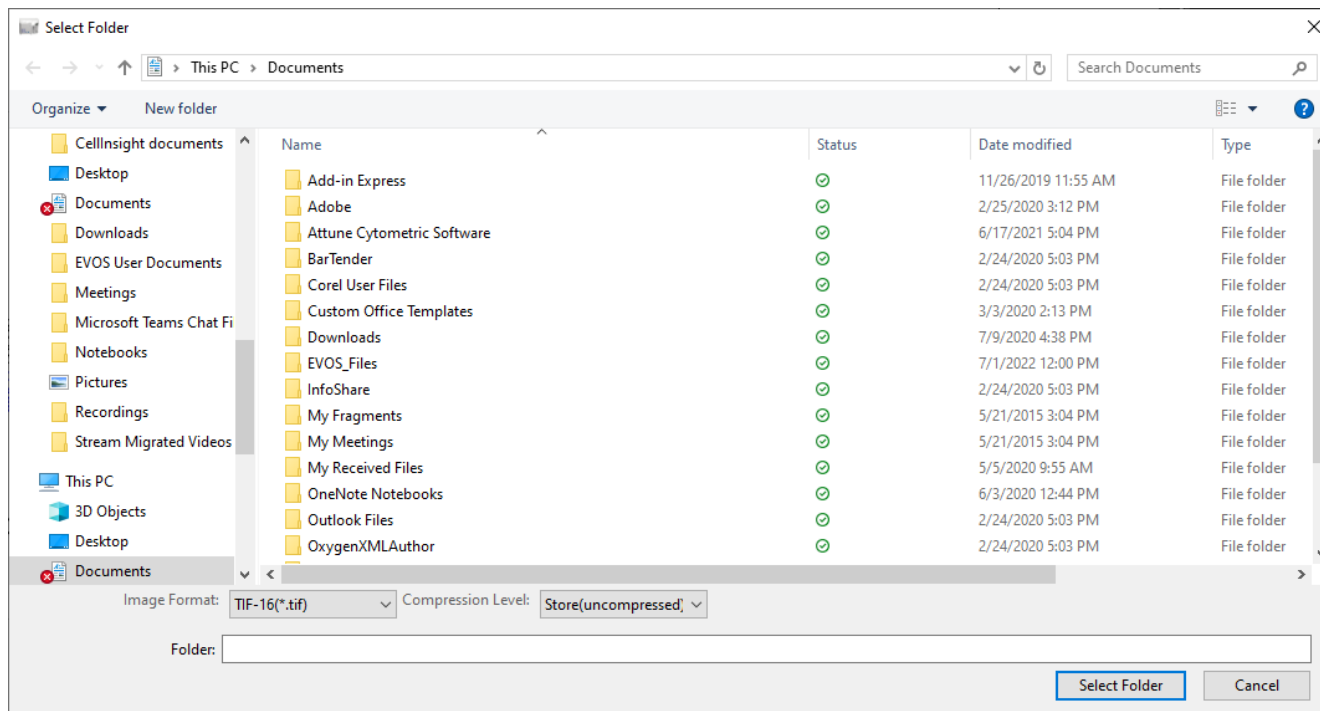
- **Export Workspace** opens the **File Save (Export)** dialog (“File Save (Export) dialog” on page 715) which enables you to select a name and location to save the Workspace file.
- The **Export Workspace** option is only enabled when a single Experiment is selected.

## Save as Template

**Save As Template** opens the Save As Template dialog (“Save As Template dialog” on page 747), which enables you to name the current Experiment and save it as a Template in the database for future reuse.

## Export Images

**Export Images** opens the **Select Folder (Folder browser) dialog** (“Folder Browser dialog” on page 723), which enables you to save the images from the selected Experiment in the desired location.



**Note:** The **Export Images** option is available only for experiments performed with an Attune™ CytPix™ Flow Cytometer, and it is active only if the experiment has image data.

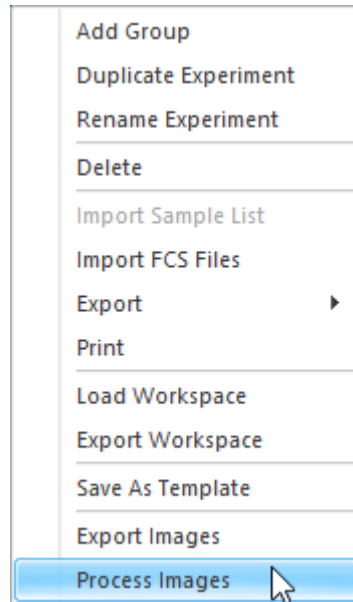
- When images are exported, all images from the selected Experiment are exported into a single zip file.
- By default, images are exported as TIF-16 files, which is the native file format collected from the Attune™ CytPix™ Flow Cytometer.
- You can select to save the images as **TIF-16**, **TIF-8**, **PNG**, **BMP**, **GIF**, **EMF**, or **JPG** files using the **Image Format** dropdown in the **Select Folder (Folder browser) dialog**.

**IMPORTANT!** When saving in any format other than **TIF-16**, images are converted to 8-bit images, which results in loss of image pixel data.

- Use the **Compression Level** dropdown to select to store the images uncompressed (**Store (uncompressed)**) or save them **Compressed** to save disk storage space.

## Process Images

**Process Images** opens the **Process Images** dialog (see Chapter 23, “Process Images dialog”), which enables you to process captured images from selected **Samples** and **Populations** using system-provided **Image Processing Models**.



- The **Process Images** option is only visible for Attune™ CytPix™ Experiments.
- You can only process images for the Experiment that is active in the Workspace. The **Process Images** option is greyed out in the **Experiment context menus** of Experiments that are not active.

---

**Note:** For more information about the Attune™ Cytometric Software image processing workflow, see the *Attune™ Cytometric Software Image Processing Workflow* (Pub. No. [MAN0028531](#)), which is available for download at [thermofisher.com](#).

---

## Group context menu

Right-click a **Group node** to show the **Group context menu**. During acquisition, all options in the **Group context menu** are disabled.

**Note:** The **Export Images** and **Process Images** options are available only for experiments performed with the Attune™ CytPix™ Flow Cytometer. These options are not visible in Attune™ NxT experiments.

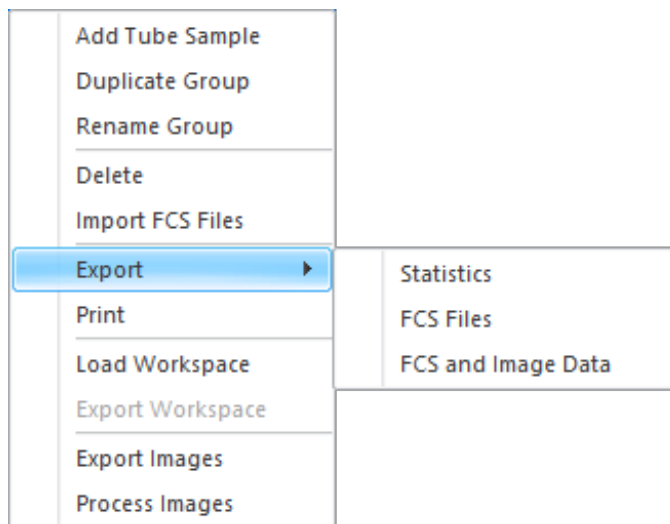


Figure 58 Group context menu

### Add Tube Sample

- **Add Tube Sample** adds a Tube sample to the selected Group.
- You can create up to 400 Samples for each Experiment. The option is disabled if there are already 400 Samples in the Experiment.
- This option is only enabled when a single Group is selected.

### Duplicate Group

- **Duplicate Group** enables you to duplicate the current Group in the current Experiment, creating an exact duplicate with the same number of Tubes and Sample-level Workspaces. If a Group includes Samples with recorded data, the duplicated Samples do not include the data.
- When a Group is duplicated, all Samples and their corresponding keyword values are duplicated. The FCS files are not duplicated.
- If duplicating a Group results in the creation of over 400 Samples for that Experiment, the new Group is created with as many Samples as possible until the 400 Sample limit is reached.
- The Duplicate option is available for Tube Experiments and for Groups in Plate Experiments that contain Tube samples. It is enabled only when a single Group is selected. Groups composed of only Well samples cannot be duplicated. If a Group includes both Well and Tube samples, the Duplicate option only duplicates the Tube samples in the Group.

- The newly created Group must have a unique name. When a Group is duplicated, it is automatically assigned a name, which consists of the same name as the current Group plus an incrementing numerical suffix. The numerical suffix starts at 1 and increments until a unique identifier is achieved.

## Rename Group

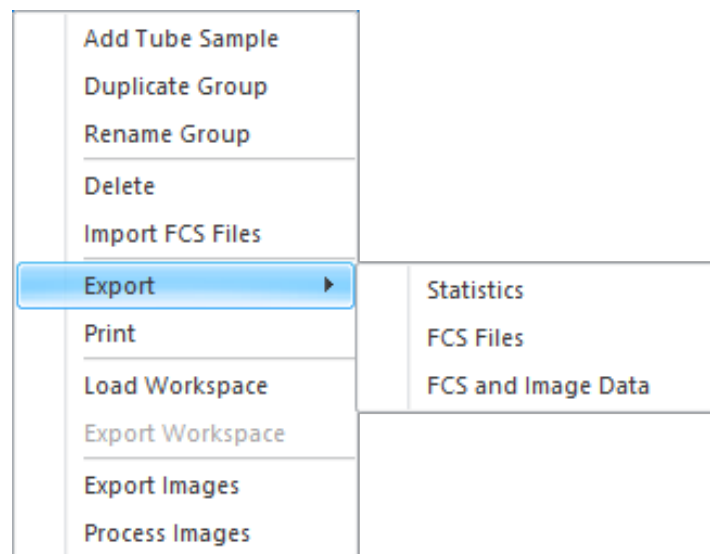
- **Rename Group** opens an edit textbox containing the name of the current Group in the edit mode.
- The **Rename Group** option is enabled when a single Group is selected.
- Clicking anywhere outside the edit textbox validates the Group name as described in “Validation of name fields” on page 301.
- You can also rename a Group by first clicking on the Group name to select it, then pressing **F2** key to enter the editing mode.

## Delete

- **Delete** opens the Deleted Items dialog (“Deleted items dialog” on page 339), which enables you to delete the selected Groups.
- The **Delete** option is only enabled when single or multiple Groups are selected in the User folder.
- If the deleted Group contains the active Sample, the software displays the Experiment-level Workspace.
- You can also use the keyboard **Delete** key to delete selected items.

## Export

**Export** displays a submenu containing **Statistics**, **FCS Files**, and **FCS and Image Data** as options.



- **Export ▶ Statistics** opens the **Export Statistics dialog** (“Export Statistics dialog” on page 737), which enables you to select Experiment-level, Group-level, or Sample-level Workspace statistics to export.
  - Statistics for all Samples in the selected Experiments are exported except Samples with no data.
  - Statistics can be exported as a single file (statistics for all Samples are combined into a single CSV file) or as individual files (statistics for all Samples are exported into separate CSV files).
  - The **Export ▶ Statistics** option is always enabled, except during acquisition.
- **Export ▶ FCS Files** opens the **File Save (Export) dialog** (“File Save (Export) dialog” on page 715), which enables you to save the selected FCS files in the desired location.
  - All FCS files in the selected Group are exported. If a Sample in the selected Group has no FCS file, then no exported FCS file is created for that Sample.
  - If the Experiment compensation settings are different from the Compensation settings contained in the FCS file, a prompt to update the compensation keyword is displayed.
  - The **Export ▶ FCS Files** option is disabled during data acquisition (Run or Record).
- **Export ▶ FCS and Image Data** opens the **File Save (Export) dialog**, which enables you to save the FCS files and imaging data together in the desired location.
  - By default, images, extended parameters (image processing data), and image mask are exported into a single zip file based on the ACS file format.
  - Exported ACS files can be imported into a Sample, which imports the FCS data and all imaging data.
  - Exported ACS files can be opened in third party software that supports the ACS standard (such as FlowJo™ Software and FCS Express™ Software).

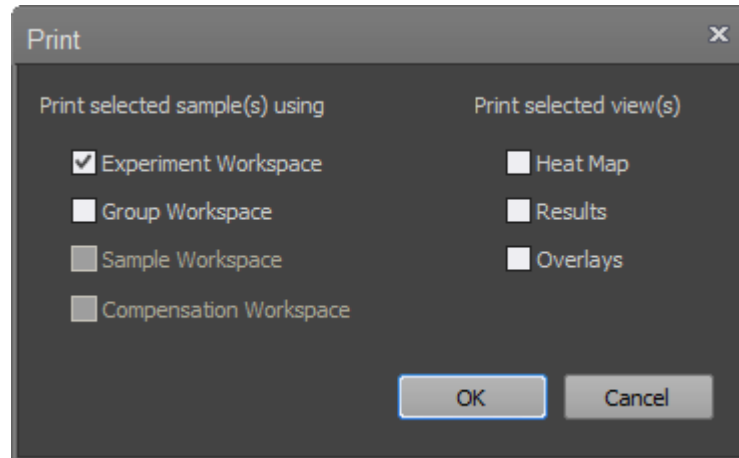
---

**IMPORTANT!** The **Export ▶ FCS and Image Data** option is available only for experiments performed with an Attune™ CytPix™ Flow Cytometer, and it is active only if the experiment has FCS and image data.

---

## Print

- **Print** opens the **Batch Print dialog** (“Batch Print dialog” on page 734), which enables you to select Workspaces (Experiment, Group, and Sample) or Views (Heat Map, Results, and Overlays) to print.
- The option to print the Compensation Workspaces is disabled, which can only be printed using the **Batch Print dialog** from the **Compensation node** (“Print” on page 337).



- The **Print** option is only enabled when single or multiple Groups in the same Experiment are selected. It is disabled during acquisition.

## Load Workspace

- **Load Workspace** opens the **File Open (Import) dialog** (“File Open (Import) dialog” on page 721), which enables you to select a Workspace file (\*.aws) to import.
- If a Workspace file is loaded, it replaces the existing Group-level Workspace.
- The **Load Workspace** option is only enabled when a single Group is selected.

## Export Workspace

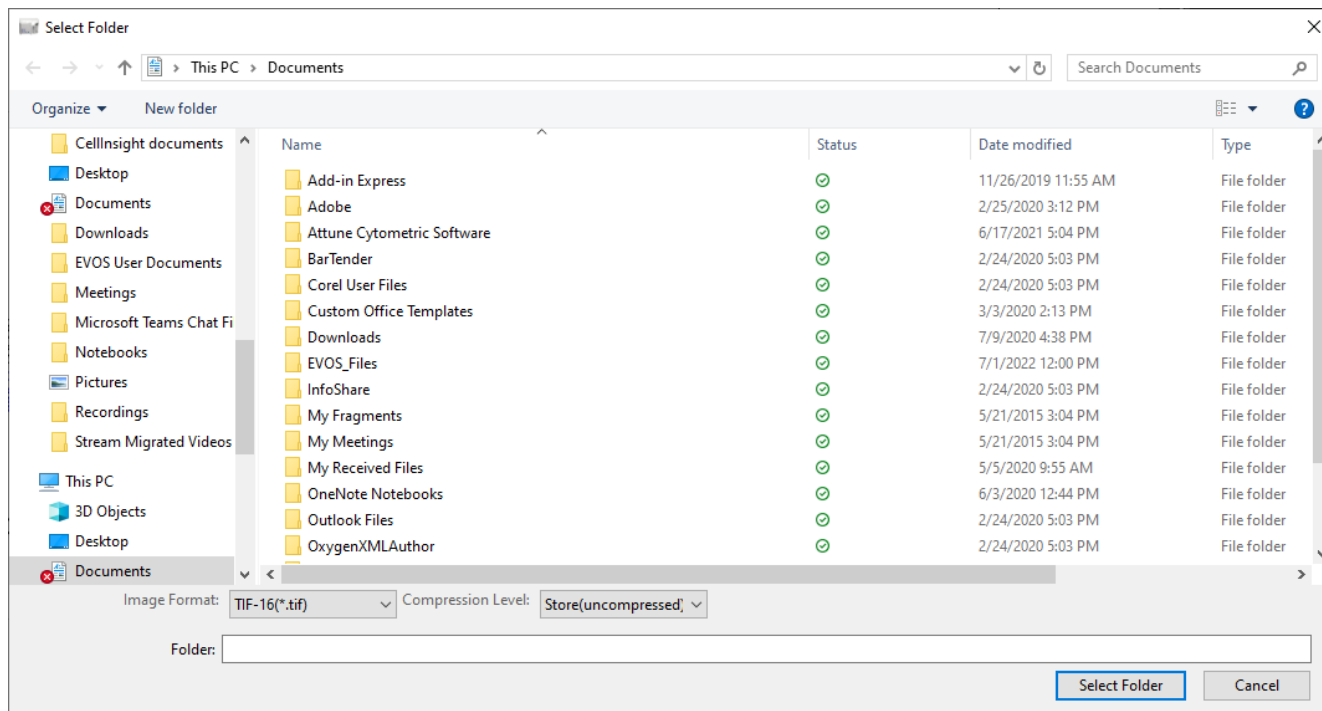
- **Export Workspace** opens the **File Save (Export) dialog** (“File Save (Export) dialog” on page 715), which enables you to select a name and location to save the Workspace file
- This option is only enabled if a Group-level Workspace exists for the selected Group.
- The **Export Workspace** option is only enabled when a single Group is selected. It is disabled when multiple Groups are selected.

## Remove Group Workspace

- **Remove Group Workspace** removes the Group-level Workspace of the selected Groups.
- This option is only available if a Group-level Workspace is present in at least one of the selected Groups.

## Export Images

**Export Images** opens the **Select Folder (Folder browser)** dialog (“Folder Browser dialog” on page 723), which enables you to save the images from the selected **Group** in the desired location.



**Note:** The **Export Images** option is available only for experiments performed with an Attune™ CytPix™ Flow Cytometer, and it is active only if the selected **Group** has image data.

- When images are exported, all images from the selected **Group** are exported into a single zip file.
- By default, images are exported as **TIF-16** files, which is the native file format collected from the Attune™ CytPix™ Flow Cytometer.
- You can select to save the images as **TIF-16**, **TIF-8**, **PNG**, **BMP**, **GIF**, **EMF**, or **JPG** files using the **Image Format** dropdown in the **Select Folder (Folder browser)** dialog.

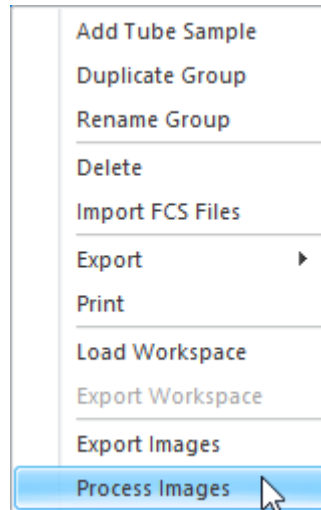
**IMPORTANT!** When saving in any format other than **TIF-16**, images are converted to 8-bit images, which results in loss of image pixel data.

- Use the **Compression Level** dropdown to select to store the images uncompressed (**Store (uncompressed)**) or save them **Compressed** to save disk storage space.



## Process Images

**Process Images** opens the **Process Images** dialog (see Chapter 23, “Process Images dialog”), which enables you to process captured images from selected **Samples** and **Populations** using system-provided **Image Processing Models**.



- The **Process Images** option is only visible for Attune™ CytPix™ Experiments.
- You can only process images for the Experiment that is active in the Workspace. The **Process Images** option is greyed out in the **Group context menus** of Experiments that are not active.

---

**Note:** For more information about the Attune™ Cytometric Software image processing workflow, see the *Attune™ Cytometric Software Image Processing Workflow* (Pub. No. [MAN0028531](#)), which is available for download at [thermofisher.com](#).

---

## Sample context menu

Right-click a **Tube sample** or **Plate sample** in **Experiment Explorer** to open the **Sample context menu**. During acquisition, all options in the **Sample context menu** are inactive.

**Note:** The **Export Images** and **Process Images** options are available only for experiments performed with the Attune™ CytPix™ Flow Cytometer. These options are not visible in Attune™ NxT experiments.

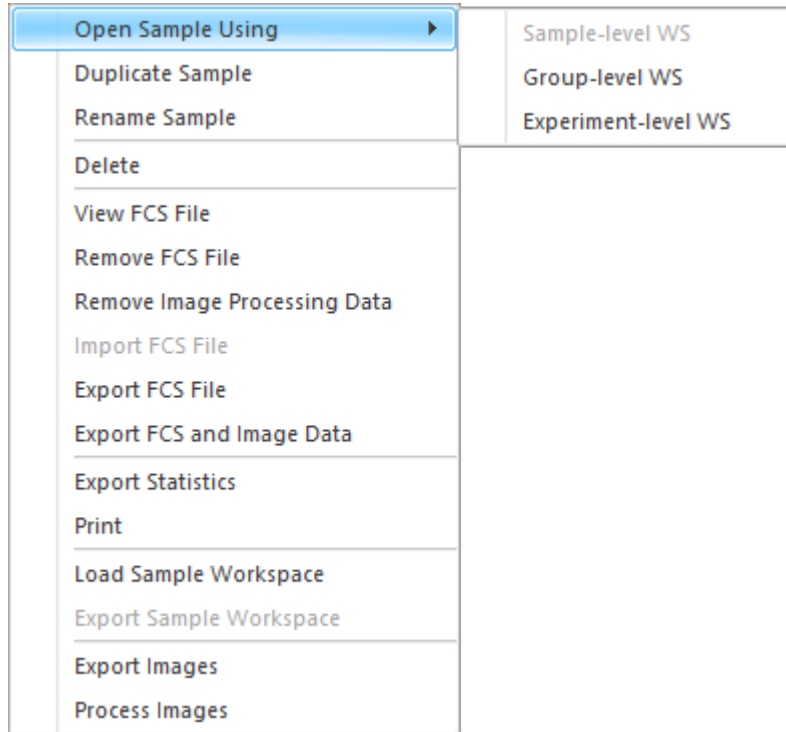
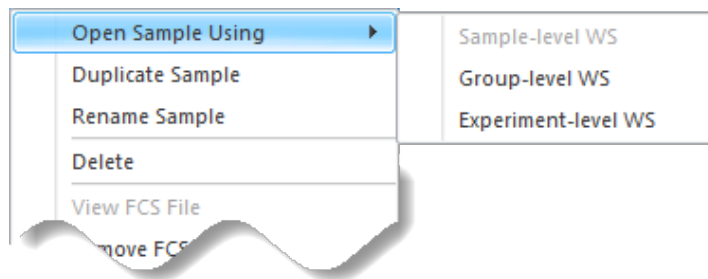


Figure 59 Sample context menu

### Open Sample Using

- **Open Sample Using** opens a secondary menu with the choice to open the selected data file using the Sample-, Group-, or the Experiment-level Workspace.



- This option enabled only when a single Sample is selected.
- The Group- and Sample-level Workspace options are enabled only if the selected Sample has a Group or Sample Workspace, respectively.

## Duplicate Sample

- **Duplicate Sample** creates a duplicate of the current Tube sample in the current Group. The Sample-level Workspace is copied, if present. All Sample-level keywords are duplicated. The Sample-level Instrument settings and the FCS file are not duplicated.
- The **Duplicate Sample** option is available only for Tube samples, and it is enabled only when a single Sample is selected.
- You can create up to 400 Samples for each Experiment. If the 400 Sample limit has been reached for the Experiment, this option is disabled.
- Each new Sample is created at the end of the current Group and must have a unique name. After a Sample is duplicated, it is automatically assigned a name, which consists of the same name as the current Sample plus an incrementing numerical suffix. The numerical suffix starts at 1 and increments until a unique identifier is achieved.
- The Sample name appears in an edit textbox ready for editing. The newly entered name is validated as described on “Validation of name fields” on page 301.

## Rename Sample

- **Rename Sample** opens an edit textbox with the name of the current Sample in edit mode.
- The **Rename Sample** option is enabled when a single Sample is selected.
- Clicking anywhere outside the edit textbox validates the Sample name as described on “Validation of name fields” on page 301.
- You can also rename a node by first clicking on the node to select it, then pressing the keyboard **F2** key to enter the editing mode.

## Delete

- **Delete** opens the **Deleted Items dialog** (“Deleted items dialog” on page 339), which enables you to delete the selected Samples.
- The **Delete** option is enabled when single or multiple Samples are selected. It is disabled if a Plate is paused or the application is in the Automation mode.
- If the active Sample is deleted, the Sample above it is activated. If there are no Samples above the deleted Sample, then the Sample below is activated. If there are no Samples, the Experiment-level Workspace is displayed. If the Experiment is deleted, the software displays the main menu.
- You can also use the keyboard **Delete** key to delete selected items.

## View FCS File

- **View FCS File** displays the FCS Sample information for the active Sample in the **FCS Sample Information panel** (Chapter 17, “FCS information panel”).
- This option is only enabled for the active Sample.

## Remove FCS File

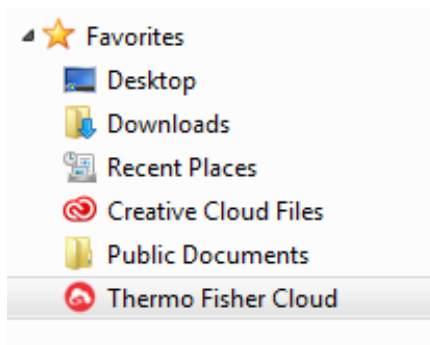
- **Remove FCS File** opens the **Deleted Items dialog** (“Deleted items dialog” on page 339).
- Confirming the **Delete** command removes the data file from the selected Samples and sends it to the recycle bin.  
Empty Sample entries are retained in the hierarchy, if the data files are removed.
- The **Remove FCS File** option is visible and enabled if one or more of the selected Samples have data present. It is disabled if a Plate is paused or the application is in the Automation mode.

## Remove Image Processing Data

- **Remove Image Processing Data** opens the **Remove Image Processing Data dialog** (“Remove Image Processing Data dialog” on page 635).
- Click **Yes** to remove the image processing data from the selected Sample and send it to the recycle bin.
- The **Remove FCS File** option is visible and active if the selected Sample has image processing data. It is inactive if a Plate is paused or the application is in the Automation mode.

## Import FCS file

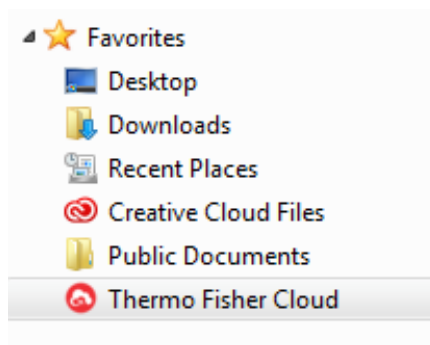
- **Import FCS File** opens the **File Open (Import) dialog** (“File Open (Import) dialog” on page 721), which enables you to select a single data file to import into the selected Sample.
- The **Import FCS File** option is visible but disabled if the selected Sample contains FCS data. It is disabled during data acquisition (Run or Record), if a Plate is paused, or if the application is in the Automation mode.
- This option is not available when multiple Samples are selected.
- If the device (instrument or software) has been registered to a Connect account, the **File Open dialogue** includes an icon for the Connect listed under **Favorites**. Clicking this icon opens a virtual folder listing all files and folders included in the Connect account. You can save the FCS file to the Connect account by selecting the desired location in the Connect account during FCS File import.



Connect account contents are only visible when viewed through the **File Open (Import)** and **File Save (Export) dialogues** from the Attune™ Cytometric Software.

## Export FCS File

- **Export FCS File** enables you to save the selected FCS file in the desired location.
- If a single Sample is selected for export, the **Export FCS File** option opens the **File Save (Export) dialog** (“File Save (Export) dialog” on page 715), which enables you to save the FCS file associated with the selected Sample using a custom name.
- If multiple Samples are selected for export, the **Export FCS File** option opens the **File Browser dialog** (“Folder Browser dialog” on page 723), which enables you to create or to select a destination folder for exporting multiple FCS files.
- If the device (instrument or software) has been registered to a Connect account, the File Open dialogue includes an icon for the Connect listed under **Favorites**. Clicking this icon opens a virtual folder listing all files and folders included in the Connect account. You can save the FCS File to the Connect account by selecting the desired location in the Connect account during Experiment export.



Connect account contents are only visible when viewed through the **File Open (Import)** and **File Save (Export) dialogues** from the Attune™ Cytometric Software.

- When exporting FCS files, you can select the FCS file name and the FCS version (FCS 3.1 or FCS 3.0). Alternatively, you can save the FCS files in the form of a CSV (.csv) file.
- If the Experiment compensation settings are different from the Compensation settings in the FCS file, the target and label names have been changed, or Experiment or custom keywords have been changed, a prompt to update the keywords is shown.

Select **Update all FCS Keywords** to update the keywords in the FCS file (for example, update the Sample name).

Select **Ignore** to close the prompt without updating the keywords.

- If a selected Sample has no FCS file, the Export FCS file option is not displayed.
- The Export FCS Files option is disabled during data acquisition (Run or Record).

## Export FCS and Image Data

**Export** ▶ **FCS and Image Data** opens the **File Save (Export) dialog**, which enables you to save the FCS files and imaging data together in the desired location.

- By default, images, extended parameters (image processing data), and image mask are exported into a single zip file based on the ACS file format.
- Exported ACS files can be imported into a Sample, which imports the FCS data and all imaging data.
- Exported ACS files can be opened in third party software that supports the ACS standard (such as FlowJo™ Software and FCS Express™ Software).

---

**IMPORTANT!** The **Export** ▶ **FCS and Image Data** option is available only for experiments performed with an Attune™ CytPix™ Flow Cytometer, and it is active only if the experiment has FCS and image data.

---

## Export Statistics

- **Export Statistics** opens the **Export Statistics dialog** (“Export Statistics dialog” on page 737), which enables you to select Experiment-level, Group-level, or Sample-level Workspace statistics to export.
- Statistics for all selected Samples that contain data are exported.
- The **Export Statistics** option is visible if one or more of the selected Samples has an associated data file. It is disabled during acquisition.

## Print

- **Print** opens the **Batch Print dialog** (“Batch Print dialog” on page 734), which enables you to select Workspaces (Experiment, Group, and Sample) or Views (Heat Map, Results, and Overlays) to print.
- The Print option is only enabled when single or multiple Samples in the same Experiment are selected. It is disabled during acquisition.

## Load Sample Workspace

- **Load Sample Workspace** opens the **File Open (Import) dialog** (“File Open (Import) dialog” on page 721), which enables you to select a Workspace file (\*.aws) to import.
- If a Workspace file is selected for import, the selected Workspace replaces the existing Sample-level Workspace.
- The **Load Workspace** option is only enabled when a single Sample is selected.

## Export Sample Workspace

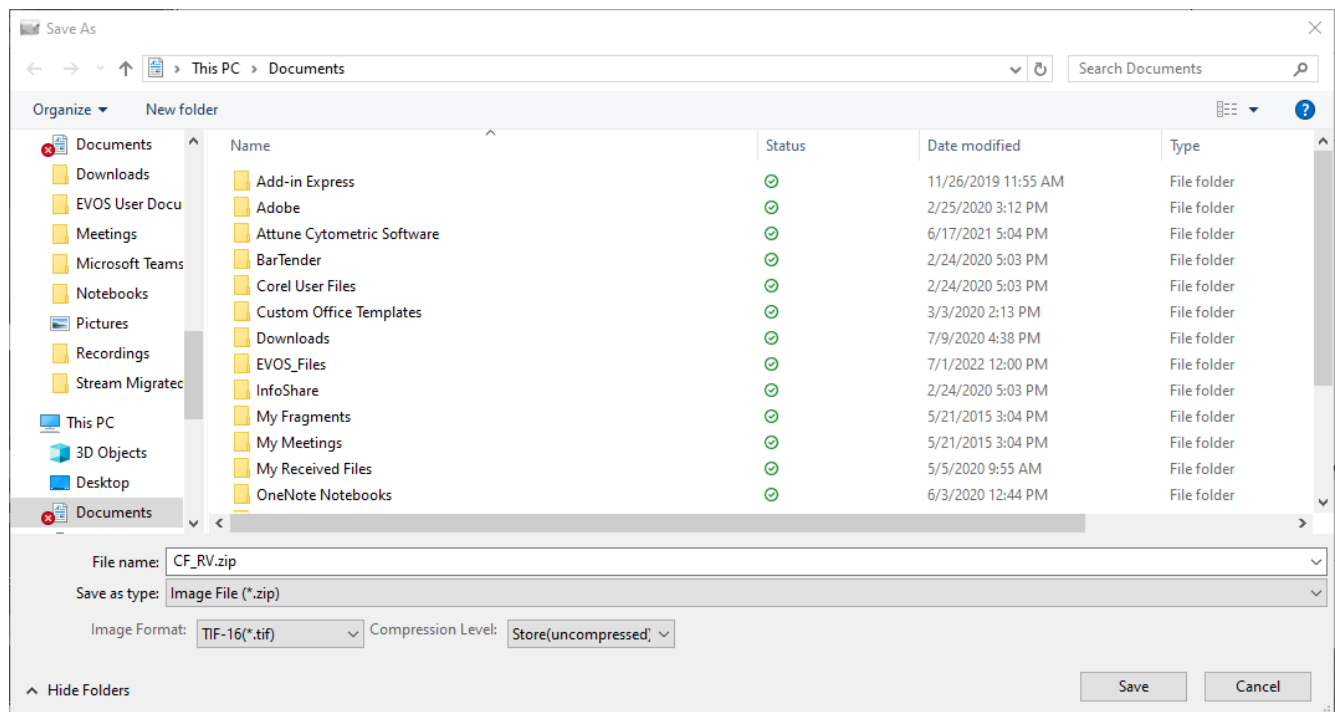
- **Export Sample Workspace** opens the **File Save (Export) dialog** (“File Save (Export) dialog” on page 715), which enables you to select a name and location to save the Workspace file.
- This option is only enabled when a single Sample is selected. It is not visible if the Sample does not contain a Sample-level Workspace.

## Remove Sample Workspace

- **Remove Sample Workspace** removes the Sample-level Workspace of the selected Samples.
- If the Sample-level Workspace that is removed is also the active Workspace, the Workspace view returns to the Experiment-level Workspace when the Sample-level Workspace is removed.
- This option is only visible if a Sample-level Workspace is present in at least one of the selected Samples.

## Export Images

**Export Images** opens the **Select Folder (Folder browser)** dialog (“Folder Browser dialog” on page 723), which enables you to save the images from the selected **Sample** in the desired location.



**Note:** The **Export Images** option is available only for experiments performed with an Attune™ CytPix™ Flow Cytometer, and it is active only if the **Sample** has image data.

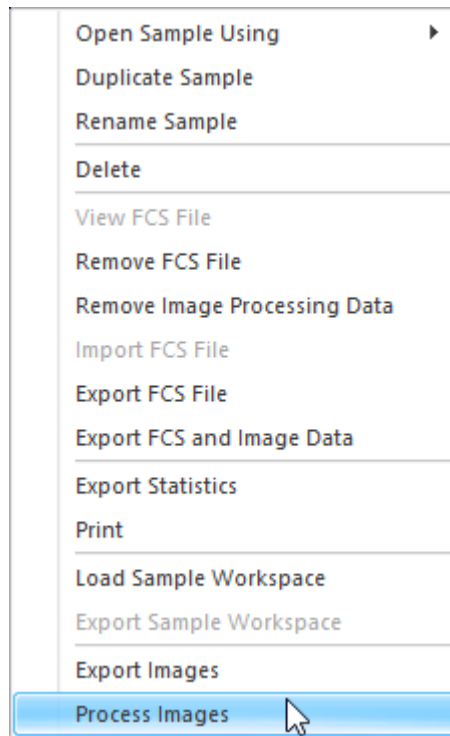
- When images are exported, all images for the selected **Samples** are exported into a single zip file.
- By default, images are exported as **TIF-16** files, which is the native file format collected from the Attune™ CytPix™ Flow Cytometer.
- Use the **Image Format** dropdown to select to save the images as **TIF-16**, **TIF-8**, **PNG**, **BMP**, **GIF**, **EMF**, or **JPG** files.

**IMPORTANT!** When saving in any format other than **TIF-16**, images are converted to 8-bit images, which results in loss of image pixel data.

- Use the **Compression Level** dropdown to select to store the images uncompressed (**Store (uncompressed)**) or save them **Compressed** to save disk storage space.

## Process Images

**Process Images** opens the **Process Images** dialog (see Chapter 23, “Process Images dialog”), which enables you to process captured images from selected **Samples** and **Populations** using system-provided **Image Processing Models**.



- The **Process Images** option is only visible for Attune™ CytPix™ Experiments.
- You can only process images for the Experiment that is active in the Workspace. The **Process Images** option is inactive in the **Group context menus** of Experiments that are not active.

---

**Note:** For more information about the Attune™ Cytometric Software image processing workflow, see the *Attune™ Cytometric Software Image Processing Workflow* (Pub. No. [MAN0028531](#)), which is available for download at [thermofisher.com](http://thermofisher.com).

---



## Compensation node context menu

Right-clicking the main **Compensation node** opens the **Compensation node context menu**. During acquisition, all options in this context menu are disabled.

**Note:** The **Export Images** and **Process Images** options are available only for experiments performed with the Attune™ CytPix™ Flow Cytometer. These options are not visible in Attune™ NxT experiments.

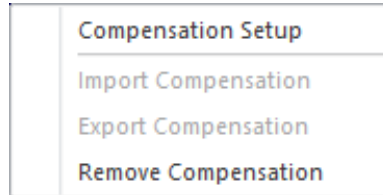


Figure 60 Compensation node context menu for Experiments without recorded Compensation (Attune™ NxT experiment)

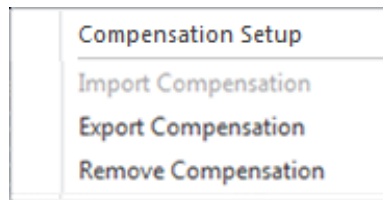


Figure 61 Compensation node context menu for Experiments with recorded Compensation (Attune™ NxT experiment)

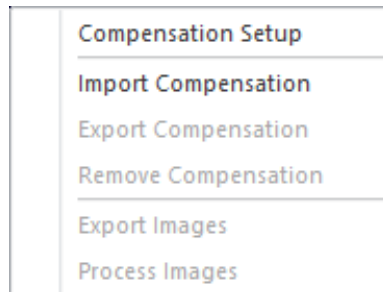


Figure 62 Compensation node context menu for Experiments without recorded Compensation (Attune™ CytPix™ experiment)

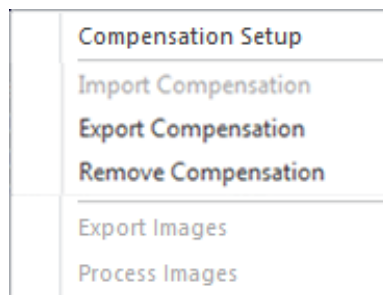


Figure 63 Compensation node context menu for Experiments with recorded Compensation (Attune™ CytPix™ experiment)

## Compensation Setup

- **Compensation Setup** opens the **Compensation Setup dialog** (“Compensation setup dialog” on page 581), which enables you to select the **Compensation Source, Compensation Measurement, Background Fluorescence Mode**, and the required compensation parameters.
- This dialog also enables you to make modifications to an existing **Compensation Setup**.
- The **Compensation Setup** option is disabled if a Plate is paused or the application is in the Automation mode. It is not visible in any Shared folder.

## Import Compensation

- **Import Compensation** opens the **File Open (Import) dialog** (“File Open (Import) dialog” on page 721), which enables you to select a Compensation Settings file (\*.acs) to import.
- Importing compensation creates Compensation controls with a read-only attribute (🔒) (“Read-only or file-based compensation” on page 332).
- The **Import Compensation** option is not enabled if any Compensation control nodes exist.

## Read-only or file-based compensation

- **Read-only or file-based compensation** is a **Compensation Setup** based entirely on the underlying Compensation XML. Read-only Compensation Controls are displayed in the Experiment Explorer with the “read-only” icon (🔒).
- A Compensation Setup cannot include both read-only controls and Sample-based controls, where an FCS file is used to determine the spillover value.
- You cannot modify read-only Compensation controls to add extra Compensation controls or to remove existing Compensation controls. You must remove the entire compensation to alter the number of controls or use Samples as Compensation controls.
- You can modify the read-only Compensation controls using any of the post-acquisition compensation modification options, including modifying the spillover values by directly typing in the Compensation matrix or using the Plot compensation modification option.
- Double-clicking on any read-only control displays the **Compensation Matrix dialog** (“Matrix dialog” on page 599).

## Export Compensation

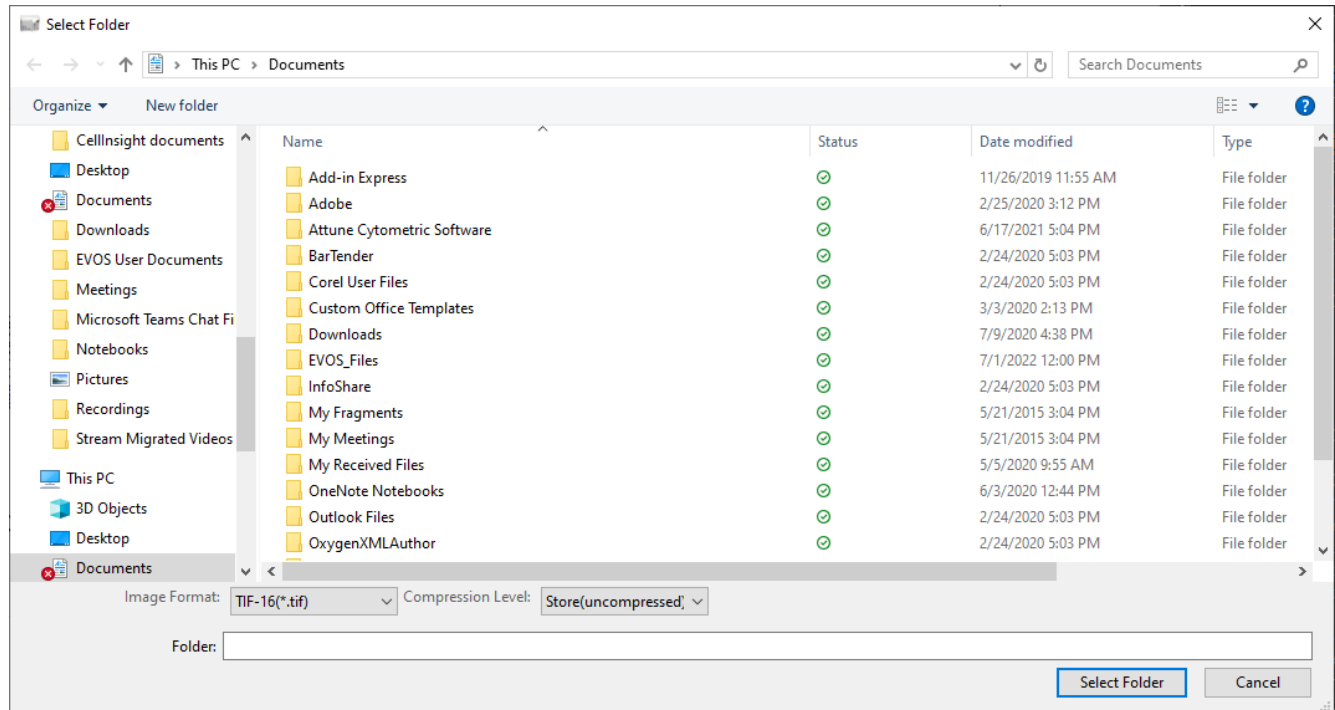
- **Export Compensation** opens the **File Save (Export) dialog** (“File Save (Export) dialog” on page 715), which enables you to select a name and location to save the Compensation Settings file (\*.acs).
- This option is enabled only when all Compensation samples have been acquired.

## Remove Compensation

- **Remove Compensation** removes the Compensation Settings and existing Compensation samples and opens the **Deleted Items dialog** (“Deleted items dialog” on page 339). If the Compensation consists of read-only controls, these are also removed.
- The **Remove Compensation** option is disabled if a Plate is paused or the application is in the Automation mode.

## Export Images

**Export Images** opens the **Select Folder (Folder browser)** dialog (“Folder Browser dialog” on page 723), which enables you to save the images from the selected **Compensation** in the desired location.



**Note:** The **Export Images** option is available only for experiments performed with an Attune™ CytPix™ Flow Cytometer, and it is active only if the **Compensation** has image data.

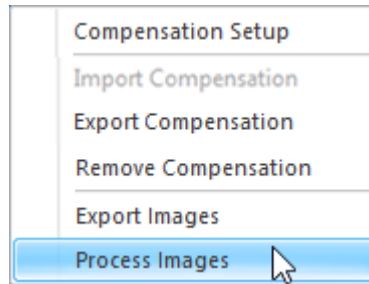
- When images are exported, all images from the selected **Compensation** are exported into a single zip file.
- By default, images are exported as **TIF-16** files, which is the native file format collected from the Attune™ CytPix™ Flow Cytometer.
- You can select to save the images as **TIF-16**, **TIF-8**, **PNG**, **BMP**, **GIF**, **EMF**, or **JPG** files using the **Image Format** dropdown in the **Select Folder (Folder browser)** dialog.

**IMPORTANT!** When saving in any format other than **TIF-16**, images are converted to 8-bit images, which results in loss of image pixel data.

- Use the **Compression Level** dropdown to select to store the images uncompressed (**Store (uncompressed)**) or save them **Compressed** to save disk storage space.

## Process Images

**Process Images** opens the **Process Images** dialog (see Chapter 23, “Process Images dialog”), which enables you to process captured images from selected **Samples** and **Populations** using system-provided **Image Processing Models**.



- The **Process Images** option is only visible for Attune™ CytPix™ Experiments.
- You can only process images for the Experiment that is active in the Workspace. The **Process Images** option is greyed out in the **Compensation node context menu** of Experiments that are not active.

---

**Note:** For more information about the Attune™ Cytometric Software image processing workflow, see the *Attune™ Cytometric Software Image Processing Workflow* (Pub. No. [MAN0028531](#)), which is available for download at [thermofisher.com](#).

---

## Compensation control context menu

Right-clicking on a **Compensation control** displays the **Compensation control context menu**. During acquisition, all options in this context menu are inactive. If the compensation is a read-only (for example, imported file), the Compensation control context menu options are all inactive.

**Note:** The **Export Images** and **Process Images** options are available only for experiments performed with the Attune™ CytPix™ Flow Cytometer. These options are not visible in Attune™ NxT experiments.

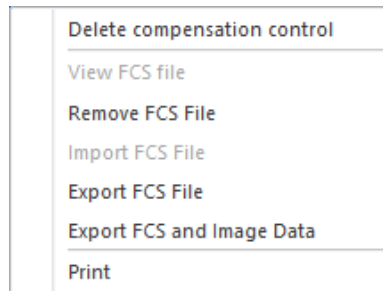


Figure 64 Compensation control context menu for controls with recorded Compensation (Attune™ NxT experiment)

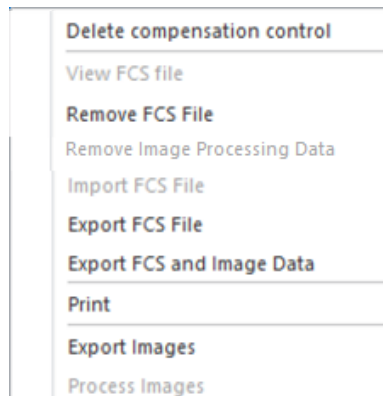


Figure 65 Compensation control context menu for controls with recorded Compensation (Attune™ CytPix™ experiment)

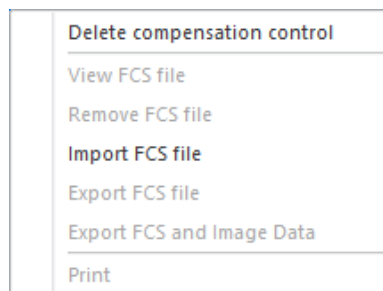


Figure 66 Compensation control context menu for controls without recorded Compensation (Attune™ NxT experiment)

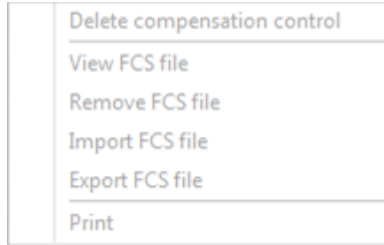


Figure 67 Compensation control context menu for read-only Compensation controls

### Delete compensation control

- **Delete compensation control** opens the **Deleted Items dialog** (“Deleted items dialog” on page 339), which enables you to delete the selected Compensation control.
- This option is disabled if a Plate is paused or if the application is in the Automation mode.
- You can also use the keyboard **Delete** key to delete selected items.

### View FCS file

- **View FCS file** opens the **FCS Sample Information panel** (Chapter 17, “FCS information panel”), which displays the FCS metadata (the text segment portion) for the selected Compensation control.
- This option is enabled only for the active Compensation control containing data.

### Remove FCS file

- **Remove FCS file** opens the **Deleted Items dialog** (“Deleted items dialog” on page 339).
- Confirming the Delete command removes the data file from the selected Compensation controls and sends the files to the recycle bin. Empty Compensation controls are retained in the hierarchy, if the files are removed.
- This option is enabled if one or more of the selected Compensation controls have data present. It is disabled, if a Plate is paused or if the application is in the Automation mode.

### Remove Image Processing Data

- **Remove Image Processing Data** opens the **Remove Image Processing Data dialog** (“Remove Image Processing Data dialog” on page 635).
- Click **Yes** to remove the image processing data from the selected Sample and send it to the recycle bin.
- The **Remove FCS File** option is visible and active if the selected Sample has image processing data. It is inactive if a Plate is paused or the application is in the Automation mode.

## Import FCS file

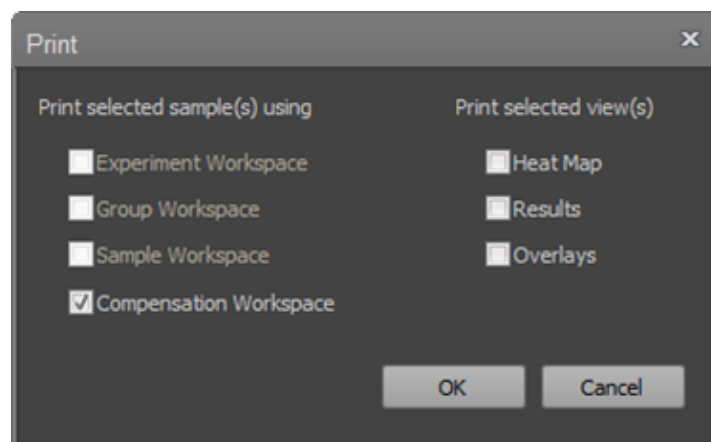
- **Import FCS file** opens the **File Open (Import) dialog** (“File Open (Import) dialog” on page 721), which enables you to select a data file to import.
- After the import of an FCS file as a Compensation control, the **Compensation Matrix** is recalculated.
- The **Import FCS file** option is visible but disabled if the selected Compensation control contains FCS data. It is disabled during data acquisition (Run or Record), if a Plate is paused, or if the application is in the Automation mode.
- This option is not visible when multiple Compensation controls are selected.

## Export FCS file

- **Export FCS file** enables you to save the selected FCS file in the desired location.
- If a single Compensation sample is selected for export, the **Export FCS file** option opens the **File Save (Export) dialog** (“File Save (Export) dialog” on page 715), which enables you to save the FCS file associated with the selected Sample using a custom name.
- If multiple Compensation samples are selected for export, this option opens the **File Browser dialog** (“Folder Browser dialog” on page 723), which enables you to create or select a destination folder for exporting multiple FCS files.  
In this case, the FCS files that are saved are named using the Experiment and the Compensation parameter name with an FCS extension.
- If a selected Sample has no FCS file, then no exported FCS file is created for that Sample.

## Print

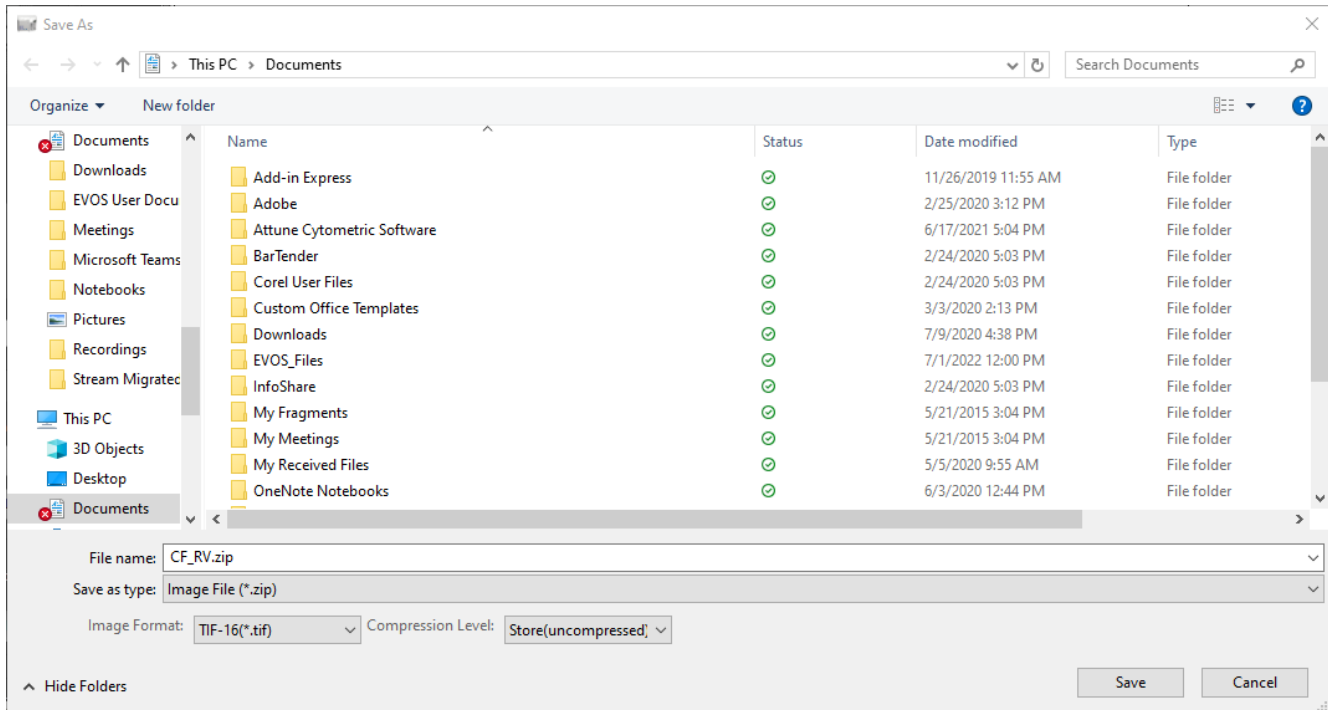
- **Print** opens the **Batch Print dialog** (“Batch Print dialog” on page 734), which enables you to select the **Compensation Workspace** and the **Heat Map, Results, and Overlays views** to print. The options to print the **Experiment, Group, or Sample Workspaces** are disabled.



- The **Print** option is only enabled when single or multiple Samples in the same Experiment compensation are selected. It is disabled during acquisition.

## Export Images

**Export Images** opens the **Select Folder (Folder browser) dialog** (“Folder Browser dialog” on page 723), which enables you to save the images from the selected **Compensation control** in the desired location.



**Note:** The **Export Images** option is available only for experiments performed with an Attune™ CytPix™ Flow Cytometer, and it is active only if the **Compensation control** has image data.

- When images are exported, all images from the selected **Compensation control** are exported into a single zip file.
- By default, images are exported as **TIF-16** files, which is the native file format collected from the Attune™ CytPix™ Flow Cytometer.
- Use the **Image Format** dropdown to select to save the images as **TIF-16**, **TIF-8**, **PNG**, **BMP**, **GIF**, **EMF**, or **JPG** files.

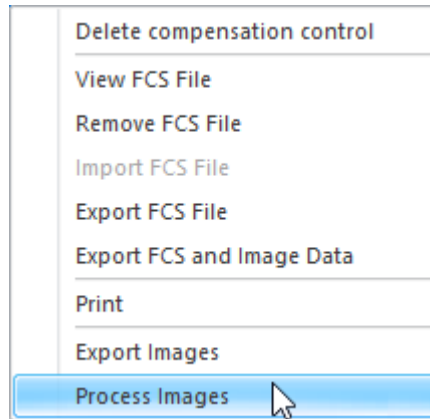
**IMPORTANT!** When saving in any format other than **TIF-16**, images are converted to 8-bit images, which results in loss of image pixel data.

- Use the **Compression Level** dropdown to select to store the images uncompressed (**Store (uncompressed)**) or save them **Compressed** to save disk storage space.



## Process Images

**Process Images** opens the **Process Images** dialog (see Chapter 23, “Process Images dialog”), which enables you to process captured images from selected **Samples** and **Populations** using system-provided **Image Processing Models**.



- The **Process Images** option is only visible for Attune™ CytPix™ Experiments.
- You can only process images for the Experiment that is active in the Workspace. The **Process Images** option is greyed out in the **Compensation control context menus** of Experiments that are not active.

---

**Note:** For more information about the Attune™ Cytometric Software image processing workflow, see the *Attune™ Cytometric Software Image Processing Workflow* (Pub. No. [MAN0028531](#)), which is available for download at [thermofisher.com](#).

---

## Deleted items dialog

The *Deleted items dialog* appears when you select **Delete** from the context menus on the Experiment Explorer or the Heat Map view, or press the **Delete key** on the keyboard to delete Tube samples, Samples, Compensation samples, Groups, Experiments, or Plates.

- All relevant prompts occur before the deletion process begins.
- Click **Yes** to move the selected FCS files to the recycle bin and remove the items from the Experiment Explorer.
- Click **Cancel** or **X** to close the dialog box (if the last item in the list) without performing an action.

### Do this for all other cases (x found) checkbox

- This checkbox appears only if multiple Experiments, multiple Groups, or multiple Samples are selected. It is unchecked by default.
- The warning displayed by the checkbox states “Do this for all other selected items (x found)”, where x is the number of selections.
- If checked, all other selected items selected are subjected to the same command.
- If left unchecked, the dialog reappears for each selected item in the queue.

## Overview

The **Collection Panel** is used for the acquisition of samples. The appearance of the **Collection Panel** varies depending on the Sample source (i.e., Plate or Tube) and the setup of the Experiment. By default, it is docked to the left of the **Main Application** area and organized into four functional groups:

- **Acquisition status** (“Acquisition status indicators” on page 344)
- **Collection** controls (“Collection controls” on page 345)
- **Collect and Display** options (“Collect and display options” on page 358)
- **Run Protocol** (“Record dialog for manual well” on page 367).

To expand or collapse the groups, click the group header. By default, only the **Run Protocol** group is expanded.

---

**Note:** Access to **Collection Panel** features can be restricted based on account permissions as described in “Options dialog – User Management options” (“User Management” on page 687) for local Attune™ Users and SAE Administrator Console (see “Roles tab” on page 886) for SAE Users.

---

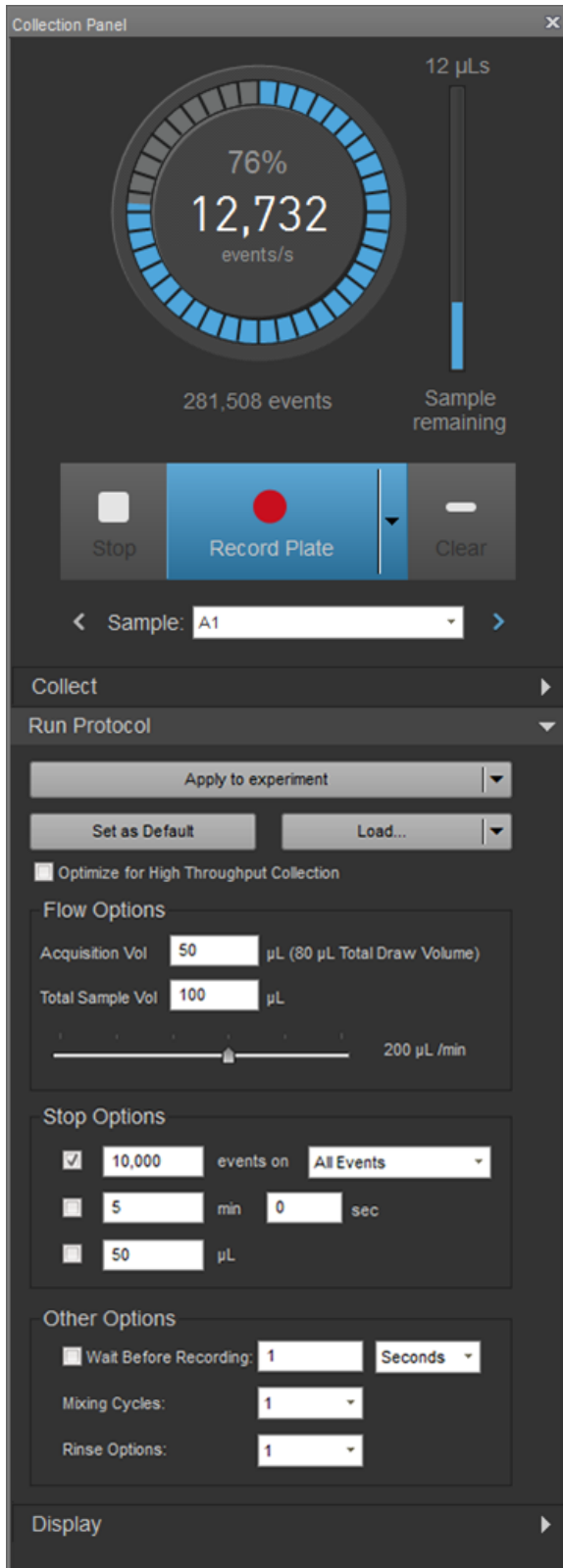


Figure 68 Run Protocol expanded (Default)

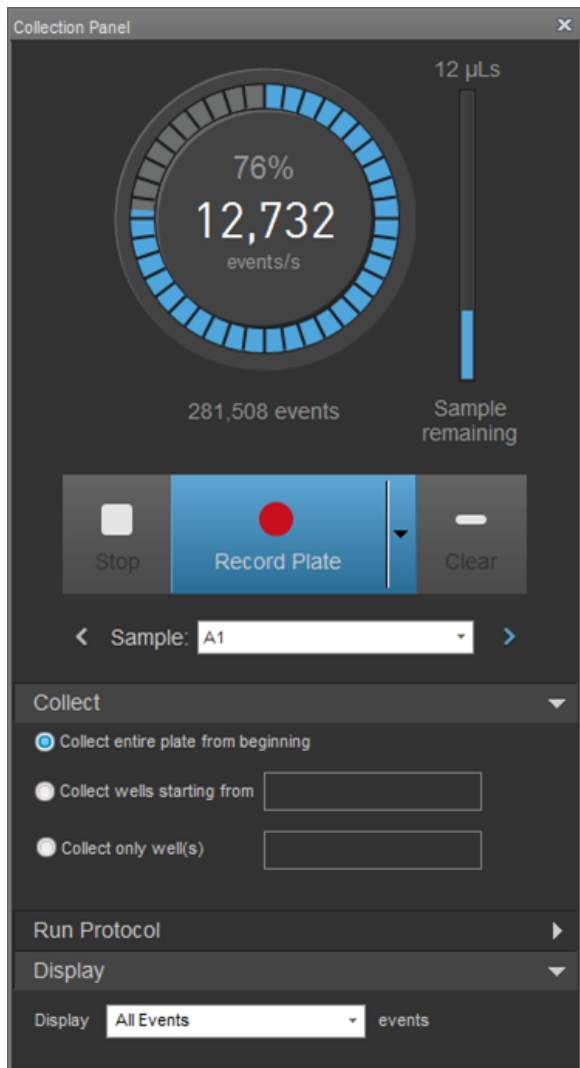


Figure 69 Collect and Display expanded

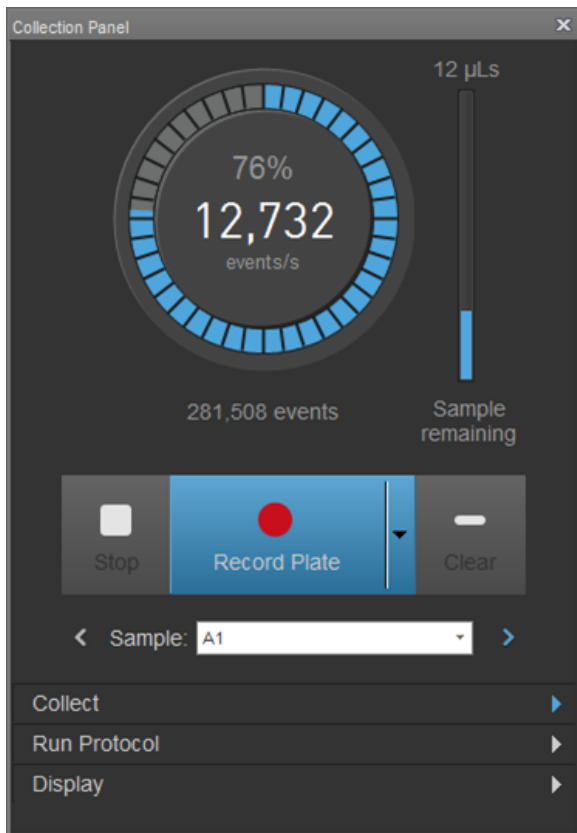
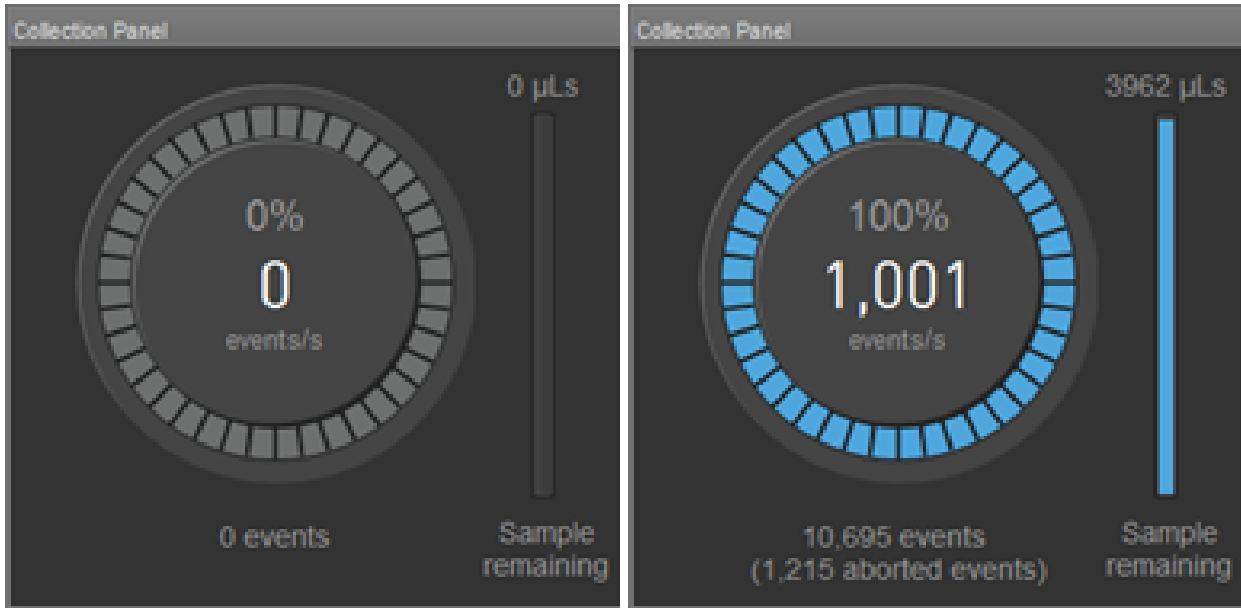


Figure 70 All collapsed

# Acquisition status

## Acquisition status indicators

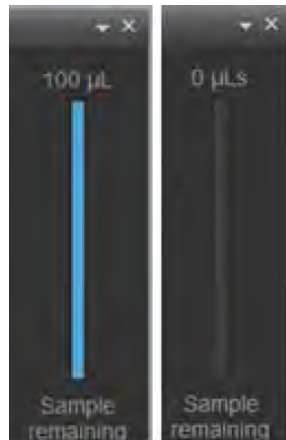
The **acquisition status indicators** provide information about the current acquisition status of the instrument.



- **Events per second (events/s)** shows the number of events currently being collected per second. This field is set to zero and does not change unless the system is acquiring data.
- **Progress dial** shows the progress of the current acquisition related to stop conditions set in the **Stop options** (“Stop Options” on page 377). The **percent progress** is displayed inside the dial above the **events per second** field.
- **Events** shows the total number of events collected. This value is reset to zero at the start of an acquisition, the start of recording the current sample, when the **Clear** button is pressed, or if a gate is moved and the stop condition is set to stop on a specific number of gated events.
- **Aborted events** shows the total number of aborted events. This value is reset to zero at the start of an acquisition, at the start of recording, or when the **Clear** button is pressed. It is only displayed if the **Exclude coincidence** checkbox is checked (“Exclude coincident events check box” on page 395).
- If **Complete Stop Condition** is selected (“Record/Save dialogs for tube sample” on page 365), the progress dial and the events field resume from where they left off.
- If multiple stop conditions are selected, the progress dial displays the condition that has progressed furthest towards being met.
- When a sample is run but not recorded, the progress dial does not increase.

## Sample remaining indicator

- The **sample remaining** indicator shows the remaining volume of the current sample.
- The **progress bar** shows the amount of sample remaining relative to the acquisition volume set as part of the Run Protocol.



## Collection controls

### Overview

The **collection controls** allow you to prompt the Attune™ instrument to run samples and record flow cytometric data.

- The **collection controls** are contextual; the options that are displayed depend on the tube or the type of well that is being acquired. Controls that are not available in a specific context are not shown or shaded gray on the panel.
- If a plate is being acquired, the default controls are the **Plate Collection Controls** (see “Collection controls – Plate” on page 350).
- If a tube is being acquired, only the **Tube Collection Controls** are available (see “Collection controls – Tube and Manual wells” on page 346).

## Collection controls – Tube and Manual wells

### Overview of collection controls for tube and manual well samples

The collection controls for Tube and Manual Well samples are similar. However, for Manual Well samples, the **Run/Record** split button has an acquisition task dropdown, and the collection controls include the **Sample navigation** and **Sample** dropdown buttons below the basic controls.

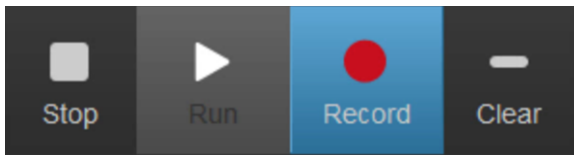


Figure 71 Collection controls for Tube samples

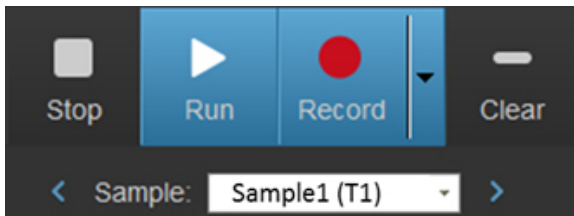


Figure 72 Collection controls for Manual Well samples

- **Run**, **Record**, **Stop**, and **Clear** buttons are the base set of buttons.
- For Manual Well samples, the **Run/Record** split button has an additional dropdown control that allows the navigation through acquisition tasks. The acquisition task dropdown is not available for Tube samples.
- When any button within the **Functions** group of the **Instrument** tab (“Functions group” on page 90) is clicked, all plate collection control buttons are inactive until the process is complete, except the navigation buttons, the acquisition task dropdown (for Manual Well samples only), and the **Save** button.
- **Run** and **Record** buttons are inactive for any current sample where the sample instrument settings do not match the system instrument configuration and for any experiment where the experiment instrument settings do not match the system instrument configuration (for example, RB vs. RBVY).



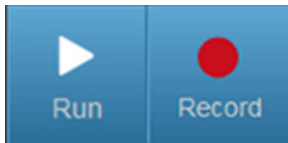
## Buttons displayed

Collection controls for Tube samples are contextual; the buttons that are displayed depend on the acquisition task and the type of Sample being acquired. The collection control buttons that are available are described below.

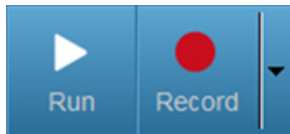
### Middle buttons

The buttons that occupy the middle section of the collection controls are used for running and recording Samples, and for saving data.

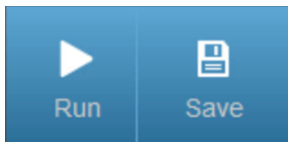
**Run/Record:** **Run** (on the left) starts the acquisition of the selected Sample; **Record** (on the right) starts the recording of the selected Sample.



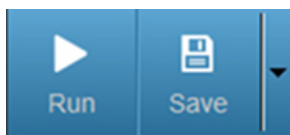
- **Run/Record** split button is displayed by default when the system is ready for acquisition.
- The **dropdown arrow** next to the center button opens the **acquisition task list**, which enables you to select an acquisition task to perform. The **acquisition task dropdown** is shown only for Manual Well samples.



**Run/Save:** **Run** reruns the selected Sample; **Save** saves the data to the FCS file for the selected Sample. If the Sample has existing data, you can append or overwrite the existing data.



- **Run/Save** split button is displayed when a Run mode acquisition has been completed for a Manual Well (that is, the end of Sample is reached or **Stop** is clicked).
- The **dropdown arrow** next to the center button opens the **acquisition task list**, which enables you to select an acquisition task to perform. The **acquisition task dropdown** is displayed only for Manual Well samples.

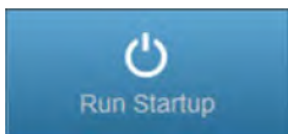


---

**Note:** The buffer size is determined by the number of events set in the **Display Events** group. By default, this is set to **All Events**.

---

**Run Startup:** Starts the **Startup** procedure.

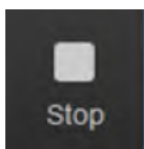


- **Run Startup** button is displayed if **Startup** has not been run before running samples or an error state requires **Startup** to be performed.

## Left buttons

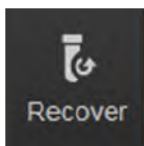
The buttons that appear on the left side of the collection controls are used for stopping sample acquisition and for recovering unused samples.

**Stop:** Stops the acquisition or recording of the current sample immediately.



- **Stop** button is available by default when the system is in acquisition state (that is, running or recording a Sample) and when running the **Startup** function.

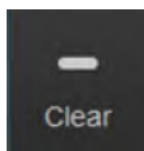
**Recover:** Recovers the unused sample to a tube.



- **Recover** button is displayed when the instrument is idle and there is a sample available to recover.
- Clicking **Recover** returns the unused sample to a tube.

## Right button

**Clear:** Clears the most recent data in the memory.



- **Clear** button is shown by default when the system is in acquisition state. It is also available post-acquisition when a sample has been overwritten in the Run mode.
- Clicking **Clear** during acquisition clears the data in memory and displays fresh data.
- Clicking **Clear** post-acquisition discards the most recent data and keeps the original data.

## Navigation buttons and sample dropdown

**Previous:** Allows navigation through previous tasks and through Samples pre and post-acquisition.



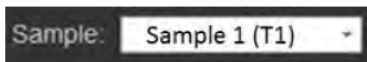
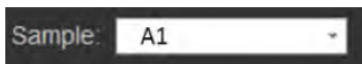
- This navigation button is available for both well and Tube samples, and can navigate from well to Tube samples, and from tube to Well samples.
- The **Previous** button is active when the current sample is not the first sample in the task list. It is inactive during acquisition.
- The button tooltip for Well samples before recording has the following form: “<phase> (sample location)”, where <phase> is Record Plate.
- After the plate is recorded, the tooltip for the wells is: “<well location>”. For example, “A1”.
- The tooltip for Tube samples has the form: “<sample name> (tube location)”, where (tube location) is T1 through T399. For example, “Sample 1 (T1)”.

**Next:** Allows navigation through samples pre and post-acquisition, and through steps in each acquisition task.



- This navigation button is available for both well and Tube samples, and can navigate from well to Tube samples, and from tube to Well samples.
- The **Next** button is active when the current sample is not the last well of the plate. However, it is active for last Tube sample and creates a new Tube sample, if clicked. The button is inactive during acquisition.
- The button tooltip for Well samples before recording has the format: “<phase> (sample location)”, where <phase> is Record Plate.
- After the plate is recorded, the tooltip for the wells is: “<well location>”. For example, “A1”.
- The tooltip for Tube samples has the form: “<sample name> (tube location)”, where (tube location) is T1 through T400. For example, “Sample 1 (T1)”.
- If a tube new sample is created by clicking **Next**, then the tooltip states: “New Sample”.

**Sample dropdown:** Allows the selection of any sample in the current acquisition task. The dropdown is inactive during acquisition.

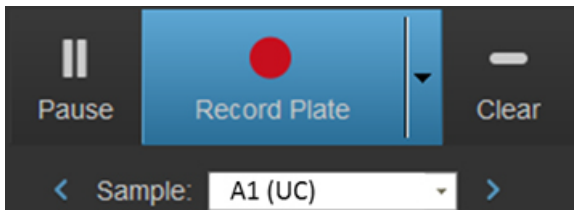


- For Sample wells, only the well location is shown. For example, “A2”.
- For Tube samples, the name is: “<sample name> (tube location)”. For example, “Sample 1 (T1)”.

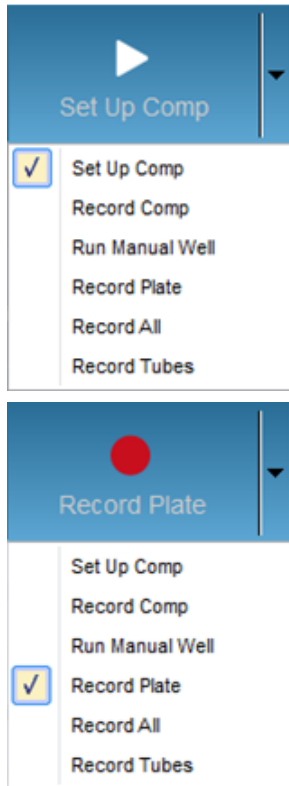
## Collection controls – Plate

### Overview of plate collection controls

**Plate collection controls** are displayed in an array of three buttons that run the basic commands for acquisition. The **Sample navigation** and **Sample** dropdown buttons are arranged below the basic controls.



- The state and the label of the buttons depend on the setup of the Plate, specifically the presence of Compensation samples, Manual Wells, and Well samples.
- The **Setup**, **Run**, and **Record** buttons are inactive for:
  - Any active Sample where the Sample-level Instrument Settings do not agree with the System Instrument Configuration.
  - Any Experiment where the Experiment-level instrument settings do not agree with the System Instrument Configuration (for example, RB vs. RBVY).
- The processing of a Plate is divided into distinct tasks. The **dropdown arrow** next to the center button opens the **acquisition task list**, which enables you to select an acquisition task to perform. The acquisition task list is available only for Plate samples.



The available tasks are:

- Set-up Compensation
- Record Compensation for the current Experiment
- Run and/or Record Manual wells for the current Experiment
- Record Plate
- Record All
- Record Tubes

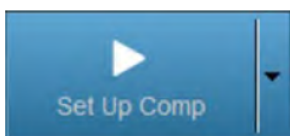
## Buttons displayed

Plate collection control buttons and their functions are described in this section. The plate collection controls are contextual; the buttons that are available depend on the acquisition task and the type of well that is being acquired. Buttons not needed in a specific context are not shown or they are inactive and shaded gray.

### Middle buttons

The buttons that occupy the middle section of the collection controls are used for running and recording samples and saving data. The labels and the visibility and functionality of the buttons depend on the plate setup and acquisition context.

**Set Up Comp:** Starts the acquisition of compensation wells in the run mode.



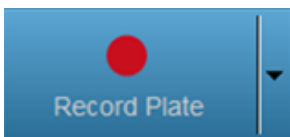
- **Set Up Comp** button is available by default if the plate contains Compensation wells and a Compensation well is active. It is also shown when **Set Up Comp** is selected from acquisition task menu.
- When the mouse pointer is hovered over the button, a tooltip states: “Allows optimization of Compensation control voltages”.

**Record Comp:** Starts recording of data from all compensation wells defined in the **Collect options** (“Collect options” on page 359).



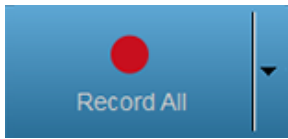
- **Record Comp** button is shown when **Record Comp** is selected from the acquisition task menu or after completing the **Set Up Comp** phase.

**Record Plate:** Starts recording of data from all Sample wells defined in the **Collect options** (“Collect options” on page 359).



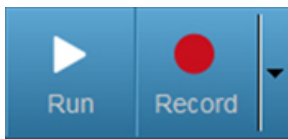
- **Record Plate** button is visible and active by default when the system is ready for recording wells and the plate contains only Well samples.
- It is also displayed when **Record Plate** is selected from the acquisition task menu or after any compensation tasks have been completed.

**Record All:** Opens the first Compensation control or Sample well in the active Plate Experiment, and automates the recording of all Compensation controls and Sample wells in the active Plate Experiment as defined in the **Collect options** (“Collect options” on page 359).



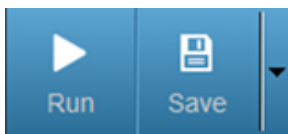
- When **Record All** is selected, you cannot manually adjust voltage settings for the Compensation controls or Manual wells.

**Run/Record:** Clicking **Run** (on the left) runs the selected Sample; clicking **Record** (on the right) records the selected Sample.



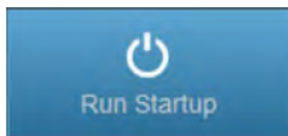
- **Run/Record** split button is displayed for Tube samples, when a Manual well is selected, when **Run Manual Well** is selected from the acquisition task menu, or after the completion of any compensation tasks (if present).

**Run/Save:** Clicking **Run** reruns the same Sample; clicking **Save** saves the data to the FCS file for the sample that has been run. If the Sample has existing data, you can append or overwrite the existing data.



- **Run/Save** split button is displayed after running a Tube sample or a Manual well.

**Run Startup:** Starts the **Startup** function.

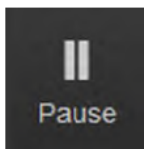


- **Run Startup** button is shown if the **Startup** function has not been run before running Samples or an error state requires **Startup** to be performed.

## Left buttons

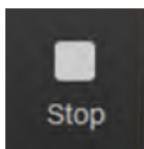
The buttons that appear on the left side of the collection controls are used for stopping and pausing sample acquisition, and for recovering preloaded samples. The visibility and functionality of the buttons depend on the acquisition context.

**Pause:** Pauses the acquisition or recording of the wells **after** the current well has finished. Pressing **Pause** during the acquisition of a well does not temporarily stop the acquisition of the well.



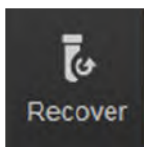
- When recording the last sample, the **Pause** button becomes a **Stop** button.

**Stop:** Stops the acquisition or recording of the wells immediately.



- **Stop** button is displayed by default during Compensation setup and when recording compensation. It is also available during **Startup**.
- When acquisition is stopped, the **Stop** button is replaced with a **Recover Sample** button if there is sample to recover.

**Recover:** Recovers the unused sample to a tube or to a well, depending on when the **Stop** button was pressed.



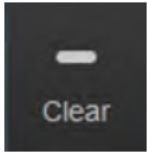
- **Recover** button is displayed when the instrument is idle and there is sample available to recover.
- Clicking **Recover** returns a preloaded sample back to a well or the unused sample to a tube.



## Right button

The right side of the collection controls only displays the **Clear** button.

**Clear:** Clears the data of the current sample and displays fresh data.



- **Clear** button is always shown. It is active when running or recording samples and wells.

## Navigation buttons

The **Previous** and **Next** navigation buttons are available for both Well and Tube samples and can navigate from Well to Tube samples and from Tube to Well samples.

**Previous:** Allows navigation through previous tasks and through Samples pre and post-acquisition.



- The **Previous** button is active when the current sample is not the first sample in the task list. The **Previous** button is inactive during acquisition.
- The tooltip for Tube samples has the form: “<sample name> (tube location)”, where (tube location) is T1 through T399. For example, “Sample 1 (T1)”.

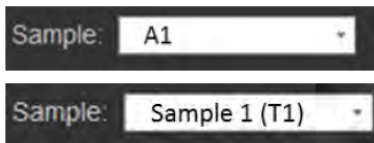
**Next:** Allows navigation through Samples pre and post-acquisition, and through steps in each acquisition task.



- The **Next** button is active when the current Sample is not the last well of the plate. However, it is active for last Tube sample and creates a new Tube sample, if clicked. The **Next** button is inactive during acquisition.
- After the Plate is recorded, the tooltip for the wells is: “<well location>”. For example, “A1”.
- The tooltip for Tube samples has the form: “<sample name> (tube location)”, where (tube location) is T1 through T400. For example, “Sample 1 (T1)”.
- If a tube new Sample is created by clicking **Next**, then the tooltip states: “New Sample”.

## Sample dropdown

**Sample** dropdown: Allows the selection of any sample in the current acquisition task. The dropdown is inactive during acquisition.



- For Sample wells, only the well location is displayed. For example, “A2”.
- For Tube samples, the name is: “<sample name> (tube location)”. For example, “Sample 1 (T1)”.

## Button availability

- The plate collection control buttons are contextual; the buttons that are displayed depend on the acquisition task and the type of well that is being acquired.
- Record Plate, Stop, and Clear buttons are the base set of buttons for plates.
- Buttons are grayed out when they are not available in a specific context.
- All plate collection control buttons are disabled if the instrument is not ready for acquisition, between wells, and after **Stop** has been clicked but the acquisition of the current sample has not been completed.
- When any button within the *Functions group* of the *Instrument tab* (“Functions group” on page 90) is clicked, all plate collection control buttons are disabled and grayed out until the process is completed, except the navigation buttons, the acquisition task dropdown, and the Save button.
- A “Bubble detected error” does not change the enabled/disabled state of a button.

## Current well indicator

- The location of the current well and the sample type that is being analyzed is displayed above the collection control buttons.
- The indicator has the format: <well location> (Compensation designation)  
For example:
  - A1 = sample in well A1
  - A1 (M) = sample in Manual well A1
  - A1 (UC) = Unstained Control sample in well A1
  - A1 (BL1) = BL1 Compensation sample in well A1

## Set up compensation workflow

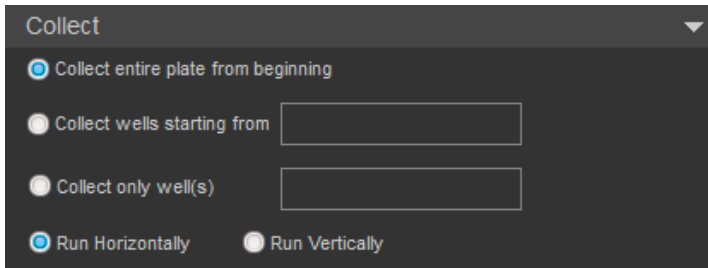
Click **Set Up Comp** to initiate the acquisition of Compensation wells.

- The order in which Compensation controls are acquired from a plate is in the same order they are listed in the Experiment Explorer:  
Unstained control, Blue laser controls, Red laser controls, Violet laser controls, and Yellow laser controls.
- Acquisition of the Compensation control continues until either **Stop** is clicked or the acquisition volume as defined by the Run Protocol (“Record dialog for manual well” on page 367) is exhausted.
- The current well remains active until the navigation buttons are used to move onto the next well within the compensation setup task or a different task is selected from the acquisition task dropdown list.
- To initiate the acquisition of the next Compensation control, you must click **Set Up Comp** again.
- Clicking the **Next** button on the last sample within the compensation setup task advances the task button to **Record Comp**.

# Collect and display options

## Collect group overview

*Collect group* allows you to control which wells are collected during acquisition. The options in this group are available only for Plate Experiments and are not displayed for Tube samples.



- Collect group contains three options for which wells to collect:
  - Collect entire plate from beginning
  - Collect wells starting from
  - Collect only well(s)

In addition, there is a selection for the Run Order, which determines the order of collection (“Run order” on page 363).

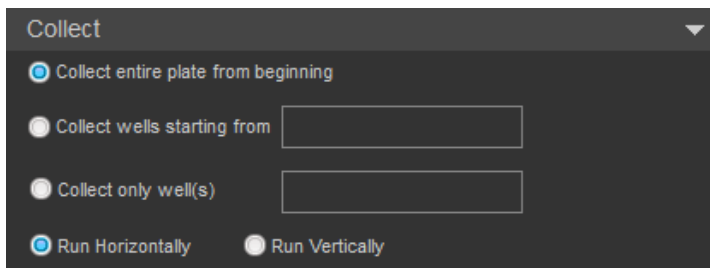
- By default, the Collect group is collapsed with the “Collect entire plate from beginning” option selected.
- The display state of Collect options persists as part of the Experiment. Collect options are also saved when duplicating, exporting, or saving a Plate as a Template, or when creating a Plate from a Template.
- If a combination of Tube and Well samples is selected in the Heat Map or in the Experiment Explorer, the options in the Collect group are disabled.
- Wells can be selected in the Heat Map to populate the “Collect wells starting from” text box.
- If the Collection panel is in the Resume state and the Collect options are modified, the Collection panel will be reset to the pre-acquisition state.
- The order of collection is based on the selected **Run Order** option (“Run order” on page 363).



## Collect options

### Collect entire plate from beginning

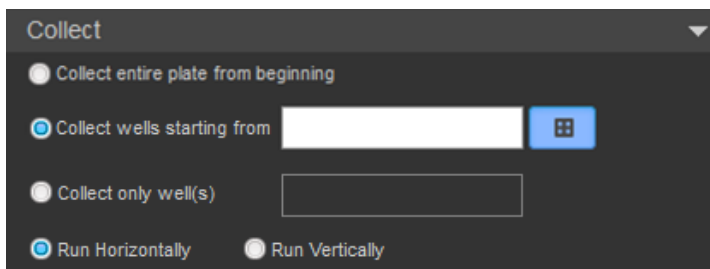
**Collect entire plate from beginning** collects all samples on a plate.



- This option is only available for Plate Experiments, and it is the default option.
- If a plate layout is specified, acquisition starts at well A1 and proceeds in the selected Run Order (“Run order” on page 363).
- The buttons that occupy the middle section of the collection controls are used for running and recording samples, and for saving data (see “Middle buttons” on page 352).

### Collect wells starting from

**Collect wells starting from** collects all samples starting at the specified well.



- To specify the starting well, enter its location in the textbox field. The entry format is the row letter followed by the column number (e.g., A1). Any letter typed is automatically capitalized.
- Alternatively, you can specify the starting well directly from the Heat Map by selecting the Heat Map interaction control (“Heat map interaction” on page 362) to the right of the textbox.



When the Heat Map interaction control is toggled on, the Heat Map of the Plate Experiment is displayed (if not already in view), allowing you to select the starting well directly from the Heat Map. When the starting well is selected on the Heat Map, the “Collect wells starting from” textbox automatically updates to display the “starting from” location.

When the Heat Map interaction control is toggled off, the selection on the Heat Map does not update the textbox contents.

- The well location entered must be a valid location and entered in the correct format. If you make an invalid entry, a warning describes the invalid entry, and the acquisition control buttons are disabled until the invalid entry is corrected. The acquisition button status indicator updates according to the well type selected.



- Invalid entries include using incorrect format (e.g., 1A), specifying a location not mapped to the selected well, entering a location not on the selected plate type (e.g., H13 on a 96-well plate), using a character that is not alphanumeric.
- Select the order of collection using the **Run Order** options (“Run order” on page 363).



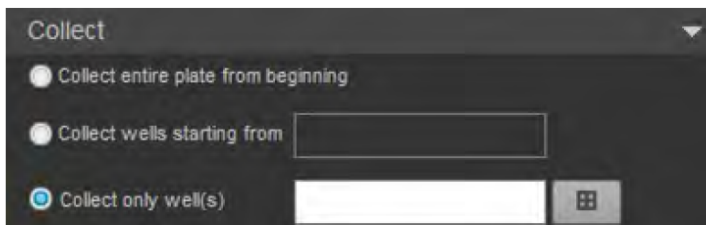
When **Run Vertically** is selected, the wells are processed from the top of the plate to the bottom of the plate, and left to right, where the collection boundary on the left is the column selected as the **Collect wells starting from**.

For example, if you select D3 as Collect all wells starting from, then E–H 1–2 are not collected, because column 3 defines the left most collection boundary.

- If a collect plate sequence is stopped, the location is automatically updated to the next well to run.

## Collect only well(s)

The **Collect only well(s)** option allows collection from the specified wells.



- To specify the wells you want to collect, enter their location in the textbox field. The entry format is one or more well locations separated by a comma or a hyphen. A well location contains a row letter followed by one or two digits for the column number (for example, B11). Any letter typed is automatically capitalized. You can use commas and hyphens for multiple entries (see below), but all other none-alphanumeric characters are invalid.
- Alternatively, you can specify the Wells you want to collect directly from the Heat Map by selecting the Heat Map interaction control (“Heat map interaction” on page 362) to the right of the textbox.



When the Heat Map interaction control is toggled on, the Heat Map of the Plate Experiment is displayed (if not already in view), allowing you to select the wells you wish to collect directly from the Heat Map as an alternative to entering the well location in the textbox.

To select multiple wells, click and drag the wells on the Heat Map or press the **Ctrl** key while clicking to select non-adjacent wells.

The Collect only well(s) textbox is automatically updated to reflect the selected wells.

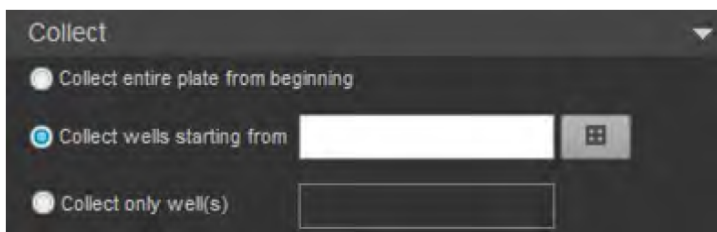
- You can enter multiple wells by using a hyphen to indicate a range of wells and a comma to separate wells and ranges. White spaces are ignored. For example: A1, B1– C12, D3.
- Where a range spans multiple rows, it is interpreted as including all wells in between. For example: B2– C6 on a 96-well plate is interpreted as containing the wells B2 to B12, C1 to C6, inclusive.
- The well location entered must be a valid location and entered in the correct format. If you make an invalid entry, a warning describes the invalid entry, and the acquisition control buttons are inactive until the invalid entry is corrected. The acquisition button status indicator updates according to the well type selected.



- Invalid entries include using incorrect format (for example, 1A), specifying a location not mapped to the selected well, entering a location not on the selected plate type (for example, H13 on a 96-well plate), using a character that is not alphanumeric.
- The order of collection is based on the order of entry of well locations. All ranges are processed left to right, then top to bottom.
- When **Collect only well(s)** is selected, the Run Order options are inactive and the wells are collected in the order they are specified in the edit control, where all ranges are processed left to right.

## Heat map interaction

When the **Collect well starting from** or the **Collect only well(s)** option is selected, a **Heat Map interaction** control becomes visible.



- When the **Heat Map interaction** button is clicked, the Heat Map is brought to the foreground.
- The textbox updates based on the wells that are selected on the Heat Map.
- For the **Collect wells starting from** option, you can select only a single well. Each time the Heat Map is clicked, the selected well is updated.
- For the **Collect only well(s)** option, you can select multiple wells by clicking and dragging to highlight the wells on the Heat Map or by pressing the **Ctrl** key while clicking to select non-adjacent wells.



## Run order

The Run Order specifies the order in which the wells are collected.

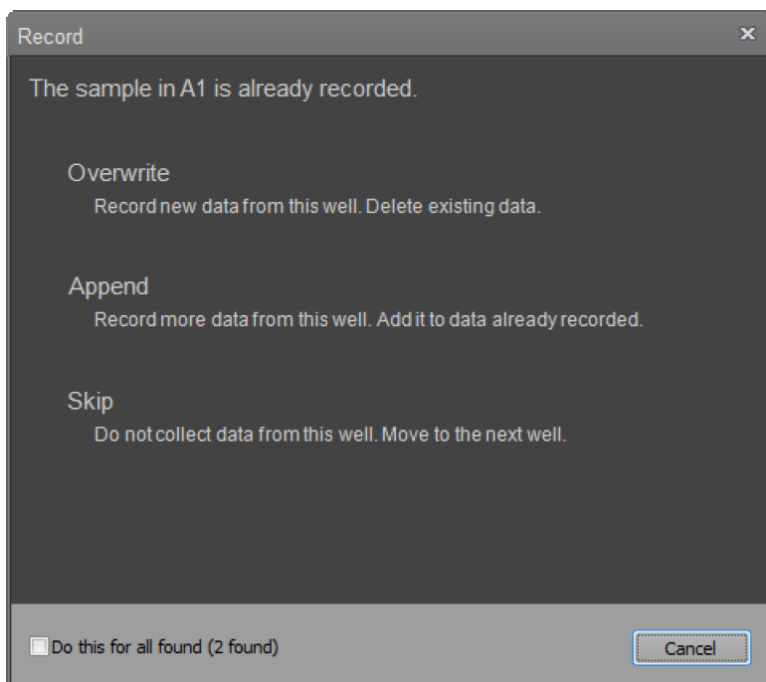


- When **Run Horizontally** is selected, wells are processed from left to right, row by row.
- When **Run Vertically** is selected, wells are processed from top to bottom, column by column.  
When **Collect wells starting from** is selected, the wells before the starting well are not collected and the well selected as the starting well defines the boundary column or row, depending on the Run Order selected.  
For example, for a plate with samples in A1–B12 with A3 as the starting well (Collect wells starting from: A3), selecting Run Horizontally results in wells A3– B12 being collected, whereas selecting Run Vertically results in wells A3–A12 and B3–B12 being collected.
- This setting only applies to the **Collect entire plate from beginning** or **Collect wells starting from** options.
- When the **Collect only well(s)** is selected, the Run Order options are disabled and the wells are collected in the order they are specified in the edit control, where all ranges are processed left to right.
- Changing the collect options does not change the selected Run Order.

## Record and save dialogs

### Record dialog for plate sample well

When **Record** is selected for a Plate sample well that has already been collected, then the **Plate Record dialog** is displayed.



The dialog heading indicates the well location and the dialog offers the following clickable options:

- **Overwrite:** Discards the data already stored in the FCS file and collects new data. Only new data are used for determining the stop condition.
- **Append:** Appends the new data to the data already stored in the FCS file. Only the new data are included in stop condition determination.  
The **Append** option is not available if the existing data file was generated in another instrument or using a different laser configuration.
- **Skip:** Skips the well without recording.
- **Do this for all found:** Applies the selection to the indicated well and all the following wells in the list that have already been collected. Otherwise the selection only applies to the well indicated in the dialog header. If there are no further wells for selection, then the next selection closes this dialog.
- **Cancel:** Closes the dialog with no action taken and the collection does not occur.

---

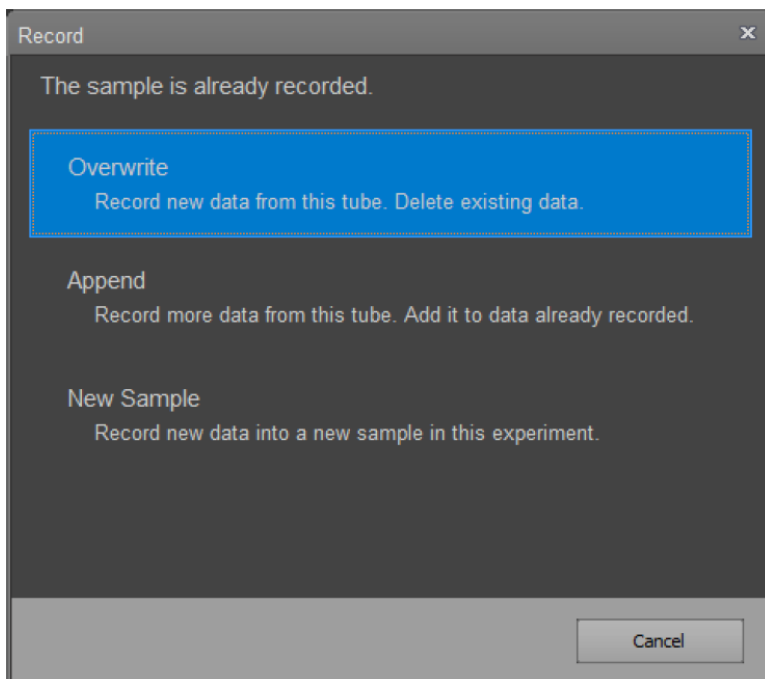
**Note:** The Run protocol for the Sample cannot be changed when overwriting, appending data, or skipping the Sample using this dialog.

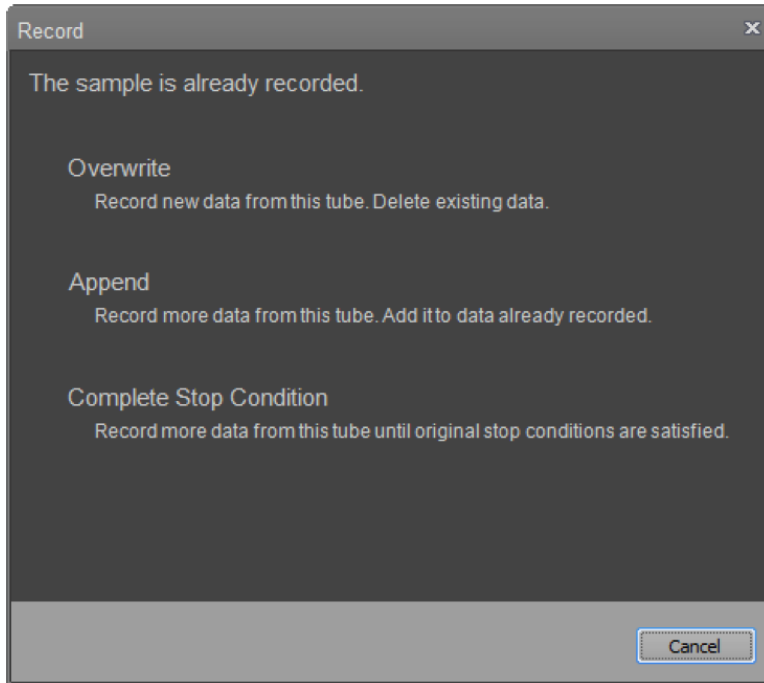
---

## Record/Save dialogs for tube sample

When **Record** or **Save** is selected for a Tube sample that already has saved data associated with it, then one of two dialogs is displayed with options to **Overwrite**, **Append**, **Complete Stop Condition**, **New Sample**, or **Cancel**.

The dialog displayed depends on the user's actions and the stop conditions. If the stop condition has already been met, the dialog on the left is displayed. If the stop condition has not yet been met, then the dialog on the right is displayed.





- **Overwrite:** Discards the data already stored in the FCS file and collects new data. Only new data are used for determining the stop condition.
- **Append:** Appends the new data to the data already stored in the FCS file. Only the new data are included in stop condition determination.  
The **Append** option is not available if the existing data file was generated in another instrument or using a different laser configuration.
- **New Sample:** Selecting this option creates a new sample in the current experiment. Data are collected for that sample using the default settings.  
The **New Sample** option is active when the **Complete stop** option is not available. It is inactive on Compensation control samples.
- **Complete Stop Condition:** Appends the new data to the data already stored in the FCS file. All data are included in stop condition determination.  
The **Complete Stop Condition** option is available only if the stop condition has not already been met, the recording was not manually stopped, and the same sample is still being run (that is, samples have not been switched).
- **Cancel:** Closes the dialog and leaves the operation in the Run mode.

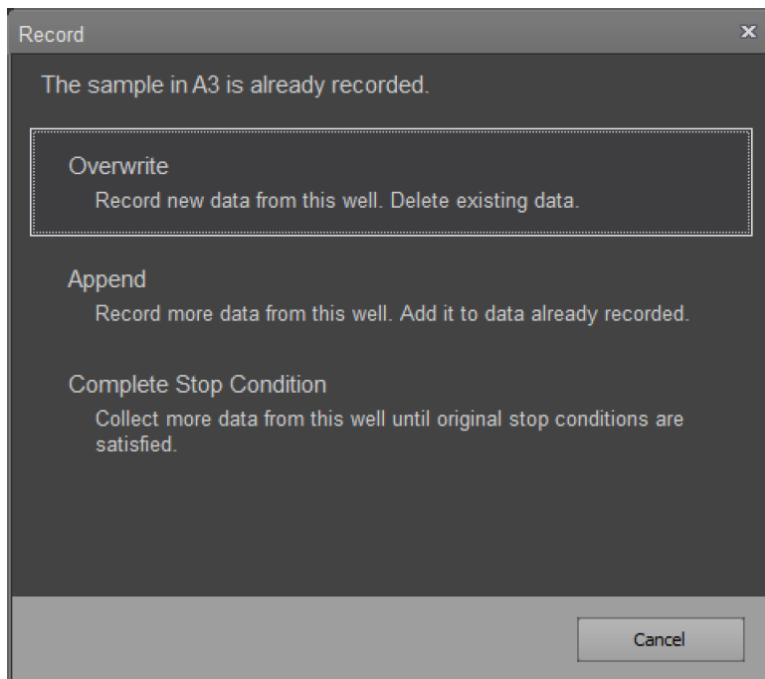
---

**Note:** The Run protocol for the Sample cannot be changed when overwriting, appending data, or skipping the Sample using this dialog.

---

## Record dialog for manual well

When **Record** is selected for a Sample well that has already been collected, then the **Plate Record dialog** is displayed.



The dialog heading indicates the well location and the dialog provides the following clickable options:

- **Overwrite:** Discards the data already stored in the FCS file and collects new data. Only new data are used for determining the stop condition.
- **Append:** Appends the new data to the data already stored in the FCS file. Only the new data are included in stop condition determination.  
This option is not available if the existing data file was generated in another instrument or using a different laser configuration.
- **Complete Stop Condition:** Appends the new data to the data already stored in the FCS file. All data are included in stop condition determination.  
This option is available only if the stop condition has not already been met, the recording was not manually stopped, and the same sample is still being run (that is, samples have not been switched).  
This option is not available if the existing data file was generated in another instrument or using a different laser configuration.
- **Cancel:** Closes the dialog with no action taken and the collection does not occur.

---

**Note:** The Run protocol for the Sample cannot be changed when overwriting, appending data, or skipping the Sample using this dialog.

---

# Run protocol

## Overview

*Run Protocol* group allows you to define the collection criteria, including the Recording and Stop options. It is also used for determining Compensation Setup options, when available. By default, the Run Protocol group is expanded.

The options displayed in this group depend on the type of sample being run and the mode of collection. The examples below show Run Protocol options for Tubes (left), Normal and Manual wells (middle), and Compensation wells (right).

**Run Protocol**

Automatically update Experiment level Run Protocol

Apply to experiment

Set as default Load...

**Flow Options**

Acquisition Vol  µL (100 µL Total Draw Volume)

12.5 25 100 **200** 500 1000 µL/min

**Stop Options**

events on

min  sec

µL

Record Events in:

Figure 73 Tube

**Run Protocol**

Automatically update Experiment level Run Protocol

Apply to experiment

Set as Default Load...

Optimize for High Throughput Collection

**Flow Options**

Acquisition Vol  µL (80 µL Total Draw Volume)

Total Sample Vol  µL

12.5 25 100 **200** 500 1000 µL/min

**Stop Options**

events on

min  sec

µL

Record Events in:

**Other Options**

Wait Before Recording:

Mixing Cycles:

Rinse Options:

Figure 74 Normal and Manual well

**Run Protocol**

Automatically update Experiment level Run Protocol

Apply to experiment

Set as Default Load...

Optimize for High Throughput Collection

**Set Up Comp Options**

Acquisition Vol  µL (80 µL Total Draw Volume)

12.5 25 100 **200** 500 1000 µL/min

**Recording Flow Options**

Acquisition Vol  µL (80 µL Total Draw Volume)

Total Sample Volume  µL

12.5 25 100 **200** 500 1000 µL/min

**Stop Options**

events on

min  sec

µL

Record Events in:

**Other Options**

Wait Before Recording

Mixing Cycles:

Rinse Options:

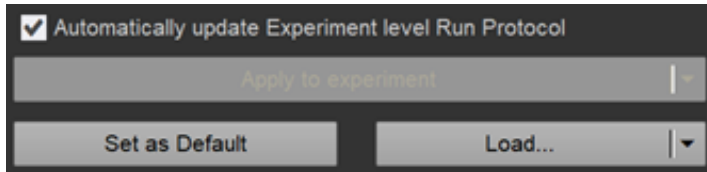
**Figure 75 Compensation well**

- Any changes made in the Run Protocol automatically apply to all active Samples, if the “Automatically update Experiment level Run Protocol” checkbox (“Automatically update experiment level run protocol” on page 371) is checked.
- If the “Automatically update Experiment level Run Protocol” checkbox is unchecked, any changes made only apply to the active Sample.
- The Run Protocols for Samples with recorded data remain unchanged. Samples with recorded data get their own Run Protocols that are not linked to the Experiment Run Protocol.



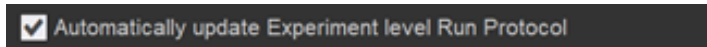
## Settings buttons

Settings buttons consist of *Apply to Experiment/Group* split button, *Set as default* button, and the *Load/Export* split button. When Run Protocol settings are modified, the changes are applied to the active Sample automatically. These buttons allow you apply the settings to other members of the current Group or Experiment.



### Automatically update experiment level run protocol

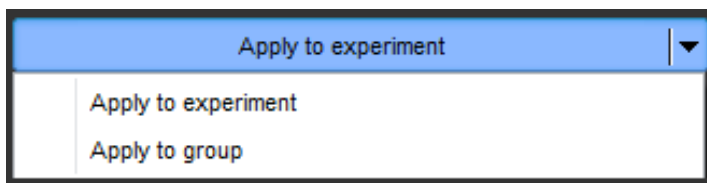
When checked, this option automatically applies the changes made in Run Protocol settings to all Samples in the Experiment that do not contain data and to all new Samples.



- By default, the “Automatically update Experiment level Run Protocol” is unchecked. Its checked/unchecked state persists as part of the user’s settings.
- If the active Sample has an associated FCS file and the Run Protocol does not match the Experiment level Run Protocol, this checkbox is disabled.
- The “Automatically update Experiment level Run Protocol” option does not apply to Compensation Controls and is disabled when a Compensation Control is active

### Apply to Experiment/Group split button

The main part of the split button contains the *Apply to Experiment* option. Clicking the dropdown arrow adjacent to the main button displays the *Apply to Group* option.



- Click **Apply to Experiment** on the main button to apply the current Run Protocol to all Samples in the current Experiment. Samples that do not have associated FCS files are updated to use the Experiment level Run Protocol.
- Select **Apply to Group** from the dropdown to apply the current Run Protocol to all Samples of the selected Group.
- When the “Automatically update Experiment level Run Protocol” is selected, this split button is disabled.
- These options are disabled if a Compensation Control is active.

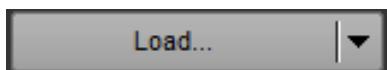
## Set as default

**Set as Default** sets the current Run Protocol to be used by default on all future Plate or Tube Experiments run by the current user.



## Load/Export split button

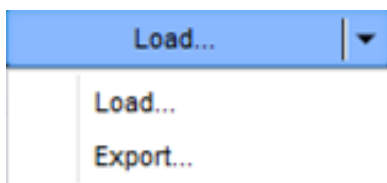
- **Load** button opens the **File Open (Import)** dialog (“File Open (Import) dialog” on page 721).



Using the **File Open (Import)** dialog, you can select a Run Protocol from a saved Run Protocol file (\*.arp).

For a sample with FCS data selecting the load option has no effect and does not open the database browser.

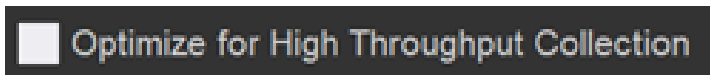
- Click the **arrow** to show the dropdown list with the **Load** and **Export** options.



- **Export** opens the **File Save (Export)** dialog (“File Save (Export) dialog” on page 715), which enables you to export and save the current Run Protocol with a user defined name as a Run Protocol file (\*.arp).

## Optimize for high throughput collection

**Optimize for high throughput collection** option optimizes the Run Protocol for high-throughput data collection.



- This option is available only when an autosampler is connected to the Attune™ instrument.
- The Optimize for High-Throughput Collection option is visible and enabled only for Plate collection (Normal well). It is visible but disabled for wells marked as Manual wells.

- When selected, the **Optimize for High-Throughput Collection** option optimizes the Run protocol for high-throughput data collection as follows:

Run protocol parameter	Attune™ NxT Auto Sampler and CytKick™ Autosampler	CytKick™ Max™ Autosampler
Stop volume	40 µL	20 µL
Acquisition volume	40 µL	20 µL
Flow rate	500 µL/minute	1000 µL/minute
Mixing cycles	1	1
Rinse between samples	1	1
Wait before recording	Unchecked	Unchecked

- If you manually change any of these settings, the checkbox is unchecked.
- If you manually adjust all settings to the conditions shown above, the checkbox becomes checked.

## Set Up Comp Options

**Set Up Comp Options** are displayed only when a Compensation well is selected. The criteria defined here are used during the set-up of Compensation wells.

### Acquisition volume

- Acquisition volume** enables you to specify the volume of sample that can be used during acquisition of a Compensation well in the Setup mode.
- The default **Acquisition volume** is 50 µL.
- The number in the **Acquisition volume** is validated on entry and must be an integer. If the number entered is outside the allowed range, it is adjusted to the nearest allowable number.
- The default, minimum and maximum values are dependent on the type of plate being used.
- The **total draw volume** (acquisition volume + dead volume) is displayed next to the acquisition volume field. This is total volume of sample drawn from the well and includes the dead volume.
- For Compensation samples, the sum of the total draw volumes for the Setup Flow options' and Recording options' acquisition volume cannot exceed the total allowable acquisition volume for the well. These limits are calculated by the software and the number entered is adjusted to the maximum allowable number.

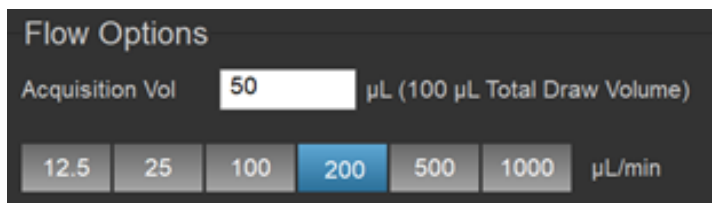
## Sample Flow Rate

The **Sample Flow Rate** controls the rate of delivery of the sample during the acquisition of Compensation samples in the Run mode.

- The flow rate values are set using the flow rate buttons. Available flow rates are:
  - 1000  $\mu\text{L}/\text{minute}$
  - 500  $\mu\text{L}/\text{minute}$
  - 200  $\mu\text{L}/\text{minute}$
  - 100  $\mu\text{L}/\text{minute}$
  - 25  $\mu\text{L}/\text{minute}$
  - 12.5  $\mu\text{L}/\text{minute}$ .
- The default flow rate is 200  $\mu\text{L}/\text{minute}$ .

## Flow/Recording Flow Options

**Flow/Recording Flow Options** are displayed for all sample types (Tube or Well samples). These settings are used when collecting data from these samples.

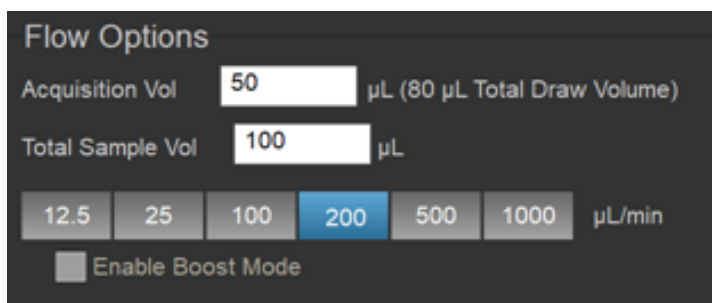


Flow Options

Acquisition Vol   $\mu\text{L}$  (100  $\mu\text{L}$  Total Draw Volume)

12.5 25 100 200 500 1000  $\mu\text{L}/\text{min}$

Figure 76 Tube Flow Options



Flow Options

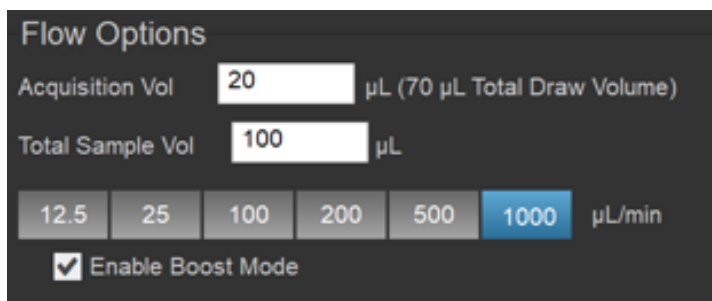
Acquisition Vol   $\mu\text{L}$  (80  $\mu\text{L}$  Total Draw Volume)

Total Sample Vol   $\mu\text{L}$

12.5 25 100 200 500 1000  $\mu\text{L}/\text{min}$

Enable Boost Mode

Figure 77 Well Flow Options



Flow Options

Acquisition Vol   $\mu\text{L}$  (70  $\mu\text{L}$  Total Draw Volume)

Total Sample Vol   $\mu\text{L}$

12.5 25 100 200 500 1000  $\mu\text{L}/\text{min}$

Enable Boost Mode

Figure 78 High-Throughput Flow Options

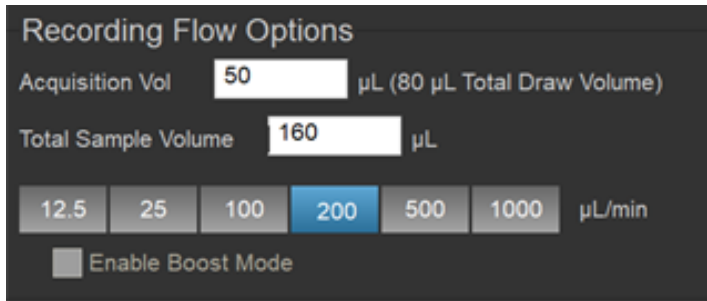


Figure 79 Recording Flow Options (for Compensation wells only)

**Note:** For Tubes and Manual wells, the group title displays **Flow Options**. For Compensation wells, the group title displays **Recording Flow Options**.

### Acquisition volume

- **Acquisition Volume** allows you to specify the volume of sample that can be used during the acquisition of the well or the tube.

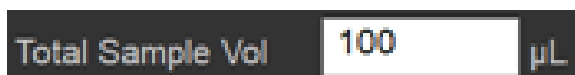


- The default Acquisition volume is 50  $\mu\text{L}$ .
- The number is validated on entry for all fields and must be an integer.
- The total draw volume (acquisition volume + dead volume) is displayed next to the acquisition volume field.
- When the acquisition volume is changed and **Automatically update Experiment level Run Protocol** or **Apply to Experiment** is selected, the changes only apply to Samples of the same type (such as wells or tube samples).

**Note:** The sample dead volume from wells is 30  $\mu\text{L}$  for sample flow rates of up to 500  $\mu\text{L}/\text{minute}$  and 50  $\mu\text{L}$  for sample flow rates of 1000  $\mu\text{L}/\text{minute}$ .

### Total Sample Volume

- **Total Sample Volume** is only displayed for Plate samples. It enables you to specify the amount of sample present in the selected well of a plate.



- The **Total Draw Volume** auto-populates the **Total Sample Volume**, but you can edit the volume to be any number greater than the **Total Draw Volume**.
- If the **Total Sample Volume** already contains a value greater than the **Total Draw Volume**, the **Total Draw Volume** does not update the **Total Sample Volume**.
- For Compensation wells, the sum of the volumes entered into the **Acquisition Volume** fields in **Setup Flow Options** and **Record Flow Options** **cannot exceed the total allowable acquisition volume** for the well. These limits are calculated by the software and the number entered is adjusted to the maximum allowable number.

---

**Note:** Setting the **Total Sample Volume** correctly ensures that Well samples are mixed adequately.

---

## Sample Flow Rate

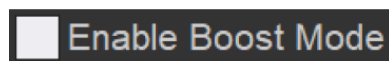
- **Sample Flow Rate** buttons control the rate of delivery of the sample during acquisition of Sample wells and Tube samples in the Record mode.



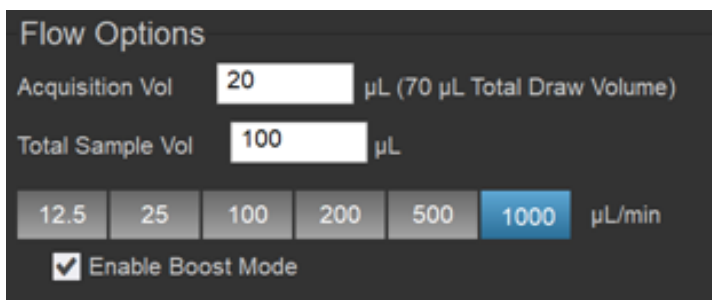
- You can select the following **Sample Flow Rates**:
  - 1000 µL/minute
  - 500 µL/minute
  - 200 µL/minute
  - 100 µL/minute
  - 25 µL/minute
  - 12.5 µL/minute.
- The default flow rate is 200 µL/minute.
- The sample dead volume from tubes is 50 µL for flow rates up to 200 µL/minute; 60 µL for 500 µL/minute, and 75 µL for 1000 µL/minute.

## Enable Boost Mode

**Enable Boost Mode** allows the processing of the Samples with reduced boost volume for high-throughput acquisition, which decreases the processing time for a standard 96-well plate from 45 minutes to 22 minutes.



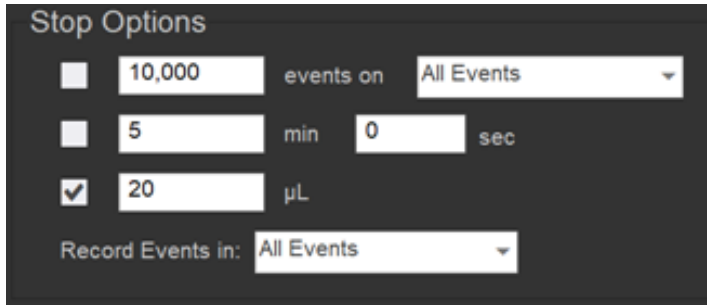
- The **Enable Boost Mode** option is available only when a CytKick™ Max Autosampler is connected to the Attune™ instrument.
- When available, you can set the **Enable Boost Mode** in the **Collection ▶ Run protocol** panel under **Flow Options**.



- The **Enable Boost Mode** option is visible only for Normal wells and Compensation wells in Plate Experiments. However, it is active only when the flow rate is set to 500 µL/minute or 1000 µL/minute.
- The **Enable Boost Mode** is not available for Manual wells or Tube samples.

## Stop Options

**Stop Options** enables you to specify when the collection of data ends. If multiple conditions are selected, the acquisition ends when any of the selected stop conditions are met. If no option is selected, the sample continues recording until the acquisition volume is exhausted.

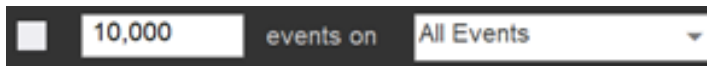


The screenshot shows the 'Stop Options' panel with the following settings:

- 10,000 events on All Events
- 5 min 0 sec
- 20  $\mu$ L
- Record Events in: All Events

### Event Count Stop

**Event Count Stop** stops acquisition when the specified number of events has been collected in the specified gate.



The close-up shows the 'Event Count Stop' control with the following settings:

- 10,000 events on All Events

- The number field enables you to enter a number between 1 and 20,000,000. The default value is 10,000. The number is validated on entry and must be an integer.
- The dropdown list displays all gates in the currently active Workspace for tubes or the Experiment Workspace for plates. An additional option, **All Events**, is always available and is the default selection.
- The following actions can result in recounting the data or changes in the **Event Count Stop**.
  - If the stop gate or a dependent gate (for example, a gate hierarchically upstream of the stop gate) is moved during acquisition, then all data are recounted. This does not occur during actual gate movement but on release of the gate to a new position.
  - If the combination of gates in the hierarchy on which the stop gate is dependent is changed during acquisition, then all data are recounted.
  - Renaming the stop gate or a dependent gate does not change the event count or the stop condition. If the stop gate is renamed, the Events dropdown is updated to show the new name when the new name is validated.
  - If the stop gate is deleted, then **All Events** option is selected and the stop option remains checked. If a dependent gate is deleted, then the stop option remains unchanged. In both cases, the data are recounted.
  - If a stop gate is undeleted as part of an undo operation, then the previous state of the **Event Count Stop** option is restored. If a stop gate or a dependent gate is undeleted, then the data are recounted.
- If recounting the data results in the stop condition being met, then the recording stops, all data are displayed, and the statistics are back-tracked to the stop event that caused the stop condition to be reached.

## Time Stop

**Time Stop** stops acquisition when the specified time has elapsed.



- You can enter a number between 0 and 59 into the **minutes** and **seconds** fields. The number is validated on entry and must be an integer. If a number greater than 59 is entered, it is adjusted to 59.
- The default stop time is 5 minutes 0 seconds.

## Volume Stop

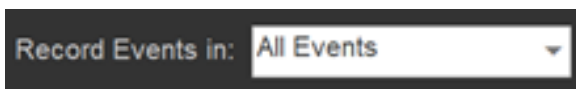
**Volume Stop** enables you to specify a delivered volume at which acquisition stops.



- The allowable range depends on the sample source, plate or tube. The number is validated on entry and must be an integer.
- If a number greater than the allowable maximum is entered, the entry is automatically adjusted. The minimum value is 1.
- For a plate, the maximum value is the entered acquisition volume.
- For a tube the maximum value is 4,000 µL.
- The default stop volume is 50 µL.

## Record Events in:

**Record Events in** allows you to set a gate where only the events that are in that gate at the end of the Sample recording are saved in the FCS file.

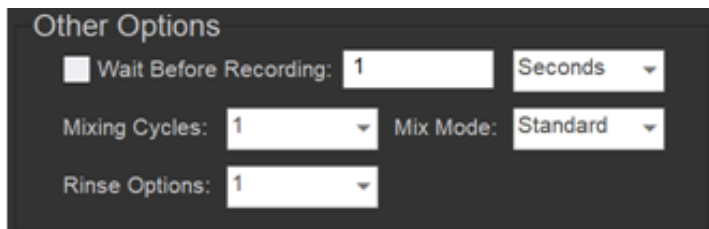


- The **storage gate dropdown** displays a list of all gates on the Experiment-level Workspace.
- By default, **All Events** is selected.
- The specified gate remains as part of the Run Protocol settings.
- As the data are acquired, **all events are displayed**, but only the events in the specified gate are saved (data exclusion only occurs on the writing of the FCS file).



## Other options

**Other Options** include **Wait Before Recording**, **Mixing Cycles**, and **Rinse Options**. These options are available when at least one well is selected.



### Wait Before Recording

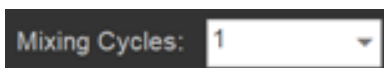
**Wait Before Recording** sets a delay for the start of acquisition.



- When the **Wait Before Recording** is enabled, the data at the start of acquisition is discarded. After the wait period or after the events count has been exceeded, data collection resumes.
- The **dropdown list** enables you to select **Seconds**, **Events**, or **Volume (µL)** to be the deciding factor for the start of recording the data.
- You can enter a number between 0 and 1,000,000 for **Events**, 0 to 30 seconds for **Time**, or 0 to 100 µL for **Volume**.
- By default, **Wait Before Recording** option is unchecked and set at 1 second.

### Mixing Cycles

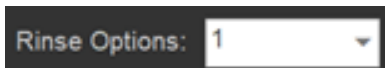
**Mixing Cycles** sets the number of mix cycles to perform on the selected wells. This option is not displayed for Tube samples.



- The **dropdown list** allows you to select between 0 and 10 mixing cycles. The default value is 1.
- The recommended number of mixing cycles is two or less to minimize the potential for creating bubbles or froth from over-mixing the samples.

## Rinse Options

**Rinse Options** sets the number of rinse cycles to perform between Well samples. This option is not displayed for Tube samples.



- The **dropdown list** enables you to select between 0 and 10 rinses. The default setting is 1.

## Mix Mode

**Mix Mode** selects between **Standard** and **Gentle** mix modes.



- The **Mix Mode** option is available only when a CytKick™ Max Autosampler is connected to the Attune™ instrument.
- When available, you can set the **Mix Mode** in the **Collection** ▶ **Run protocol** panel under **Other Options**.



- The **Mix Mode** option is available only when a Plate well is selected. It is not available for Tube samples.
- When the option is set as **Standard**, the Sample is mixed using the normal mixing speed.
- When the option is set to **Gentle**, the Sample is mixed and aspirated at a slower speed.

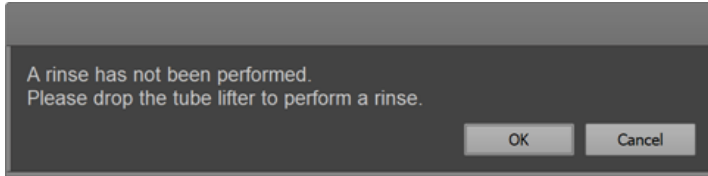
---

**Note:** For fragile cells, viscous samples, or samples prepared in viscous buffers, use the **Gentle** mix mode.

---

## Rinse reminder dialog

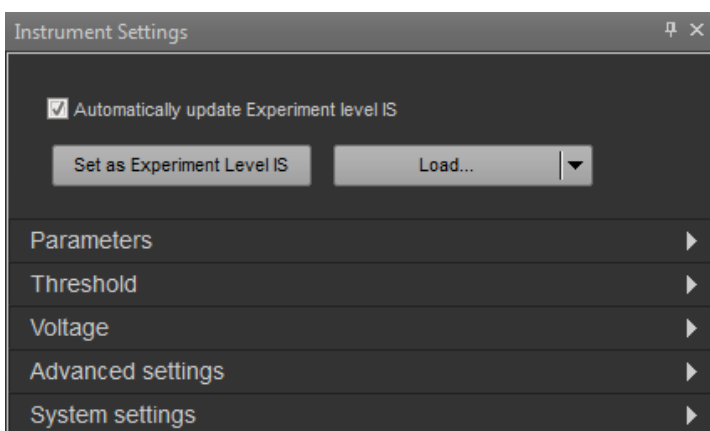
**Rinse reminder dialog** is displayed if you have acquired a sample from a tube and have not dropped the tube lifter to start an automatic rinse within 3 minutes of sample acquisition.



- Click **OK** to close the dialog. The dialog is presented again after another 3 minutes if the tube lifter was not dropped.
- Drop the tube lifter to automatically close the dialog and perform an automatic rinse.
- If you run another acquisition from the same tube within the 3 minutes, the timer is stopped and reset at the end of acquisition.

## Overview

*Instrument Settings* panel allows you to edit acquisition parameters, create custom parameters, define threshold and voltage settings, and edit advanced instrument settings. By default, it is docked to the left of the Main Application area (“Main application workspace” on page 56).



The Instrument Settings panel is organized into six functional groups in the following order:

- *Default and Load options* (“Default and load options” on page 383)
- *Parameters* (“Parameters” on page 385)
- *Threshold* (“Threshold” on page 392)
- *Voltage* (“Voltage” on page 396)
- *Advanced™ settings* (“Advanced™ settings” on page 398)
- *System settings* (“System instrument settings” on page 401)

The Instrument Settings options are contextual; controls that are not available in a specific context are not shown or shaded gray on the panel.

## General panel properties

- The docking properties of the Instrument Settings panel are described on “Docking locations” on page 58.
- If the instrument settings for the active sample do not match the system instrument configuration, the instrument settings panel displays the following message:  
“Instrument settings cannot be displayed for the selected sample. The instrument configuration is not supported”.
- The same warning message is displayed for any Experiment where the instrument settings for the Experiment do not match the system instrument configuration (e.g., RB vs. RBVY). Contents of other instrument settings panels are not displayed.
- Typing into the textbox and pressing **Enter** updates the value in the textbox. Pressing **Tab** updates and moves to the next item specified.

## Default and load options

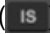
### Overview

*Default and Load options* consist of the three controls that lie at the top of the Instrument Settings panel. They allow automatic updating of Experiment-level instrument settings, one-time updating of Experiment-level instrument settings, and saving of instrument settings as default.

### Automatic update option

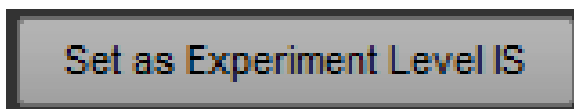
**Automatically update experiment instrument settings** checkbox enables automatic update of Experiment-level instrument settings to the previously recorded Sample-level instrument settings for each Sample when instrument settings are changed.


- The checkbox works for Tube samples only, and it is checked by default.
- Compensation samples (Tube or Well) always update the Experiment-level instrument settings. When Compensation samples are active, the checkbox is automatically checked and inactive (only when running compensation).
- Behavior, if the checkbox is checked:
  - If the instrument settings are modified, the Experiment-level instrument settings are automatically updated.
  - If the instrument settings are modified on subsequent Samples, the Experiment-level instrument settings are automatically updated.
  - If any Sample has recorded data, change of the Experiment-level instrument settings result in the recorded Samples displaying the **Instrument Settings badge** (IS), which indicates that the Sample now has Sample-level instrument settings independent of the Experiment. The Sample-level instrument settings reflect the instrument settings used at the time the Sample was recorded
  - Changes to Instrument Settings after a Sample has been recorded that alters the state of the **IS** badge is a change to the Parameter label (target and label) after the Sample is recorded.

- Behavior, if the checkbox is unchecked:
  - If the active Sample's instrument settings are modified, the modified settings do not update the Experiment-level IS. Any additional changes to the instrument settings only apply to the active Sample.
  - The Sample displays the **IS** badge (  ), which indicates that a Sample-level instrument settings has been created independent of the Experiment-level instrument settings.
  - When the checkbox is unchecked, any change of the instrument settings of the current Sample only affect the current Sample.
  - If a Sample-level instrument settings is created for a Sample, the checkbox is unchecked and inactive until the Sample's settings are applied to the Experiment either by using the **Set as Experiment Level IS** button or by dragging the **IS** badge to the Experiment level.

## Set as experiment IS

- Clicking **Set as Experiment Level IS** for Tubes saves the Instrument Settings of the current active Sample as the Experiment-level Instrument Settings.



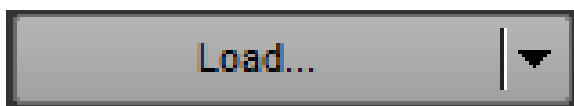
- Experiment-level Instrument Settings are indicated by the **IS** badge (  ), which is displayed to the right of the Experiment name in Experiment Explorer.
- Subsequent Samples in this Experiment will have these Instrument Settings as default.
- If the Experiment contains Samples with recorded data, those Samples show the **IS** badge

## Save as Default split button

- The **Save as Default** button allows the Instrument Settings to be saved as a default.



- The split button is inactive during tube acquisition, well acquisition, or when processing a plate.
- **Save as Default:** Saves all changes to the Instrument Settings as default Instrument Settings to be used for all new Experiments. This includes all active and selected parameters and measurements, targets, fluorophores, thresholds, voltages, custom parameters, and advanced settings.
- **Load:** Opens the **File browser** dialog ("Folder Browser dialog" on page 723).



Using the **File browser** dialog, you can select an Instrument Settings file (\*.ais) from a saved location

The **Load** button is available only when the Experiment does not include data.

- **Export:** Opens the **File Save (Export)** dialog (“File Save (Export) dialog” on page 715). Using the **File Save (Export)** dialog, you can save the current Instrument Settings with a user-defined name in the specified folder.  
You will be prompted for confirmation before overwriting an existing Instrument Settings file.

## Parameters

### Overview

**Parameters** group enables you to select the parameters to be recorded in the Experiment, based on the selected instrument configuration.

- Parameters group contains the following controls:
  - **Configuration** dropdown (“Configuration dropdown” on page 387)
  - **Parameter selector** checkboxes (“Parameter selector” on page 389)
  - **Target** textbox (“Target textbox” on page 390)
  - **Fluorophore** textbox (“Label textbox” on page 391)
- Parameters are listed in the following order:
  - Scatter channels (FSC, SSC)
  - Blue laser channels (BL)
  - Red laser channels (RL)
  - Violet laser channels (VL)
  - Yellow laser channels (YL)

Parameters

Baseline/PT Config BRVY

Time Update all samples

Event

Enabled	Target	Label	A	H	W
<input checked="" type="checkbox"/>			<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	FSC		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	SSC		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	BL1		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	BL2		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	BL3		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	RL1		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	RL2		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	RL3		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	VL1		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	VL2		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	VL3		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	VL4		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	YL1		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	YL2		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	YL3		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	YL4		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>



- Only lasers and channels in the installed laser configuration are available in the Parameters options. The list for possible configurations and their respective laser and channel order is shown in the following table.

B	BR	BY	BG	BV	BV6	BRY	BRG	BVY	BVG	BRV	BRV6	BRVY	BRVG	BRV6Y
FSC of B	FSC of B	FSC of B	FSC of B	FSC of B	FSC of B	FSC of B	FSC of B	FSC of B	FSC of B	FSC of B	FSC of B	FSC of B	FSC of B	FSC of B
SSC of B	SSC of B	SSC of B	SSC of B	SSC of B	SSC of B	SSC of B	SSC of B	SSC of B	SSC of B	SSC of B	SSC of B	SSC of B	SSC of B	SSC of B
BL1	BL1	BL1	BL1	BL1	BL1	BL1	BL1	BL1	BL1	BL1	BL1	BL1	BL1	BL1
BL2	BL2	BL2	BL2	BL2	BL2	BL2	BL2	BL2	BL2	BL2	BL2	BL2	BL2	BL2
BL3	BL3	BL3	BL3	BL3	BL3	BL3	BL3	BL3	BL3	BL3	BL3	BL3	BL3	RL1
BL4	BL4	YL1	GL1	BL4	VL1	RL1	RL1	VL1	VL1	BL4	RL1	RL1	RL1	RL2
	RL1	YL2	GL2	VL1	VL2	RL2	RL2	VL2	VL2	RL1	RL2	RL2	RL2	RL3
	RL2	YL3	GL3	VL2	VL3	RL3	RL3	VL3	VL3	RL2	RL3	RL3	RL3	VL1
	RL3	YL4	GL4	VL3	VL4	YL1	GL1	VL4	VL4	RL3	VL1	VL1	VL1	VL2
				VL4	VL5	YL2	GL2	YL1	GL1	VL1	VL2	VL2	VL2	VL3
					VL6	YL3	GL3	YL2	GL2	VL2	VL3	VL3	VL3	VL4
						YL4	GL4	YL3	GL3	VL3	VL4	VL4	VL4	VL5
								YL4	GL4	VL4	VL5	YL1	GL1	VL6
											VL6	YL2	GL2	YL1
												YL3	GL3	YL2
												YL4	GL4	YL3

- If you change the instrument configuration, target, or fluorophore for any sample post-recording, a **notification badge** (🔔) is displayed next to the affected parameter or the **Configuration** dropdown (“Notification badge” on page 391), and a notification key becomes visible at the bottom of the panel.
- Clicking the **notification badge** reverts the change to the value stored in the FCS file.

## Configuration dropdown

**Configuration** dropdown enables you to select a filter configuration and update the choice of fluorophore for each enabled parameter for the selected configuration.



- The **Configuration** dropdown list is linked to the filter configuration manager and contains all available configurations saved by the current user.
- The configuration selected from the dropdown list updates the choice of fluorophores for each enabled parameter. The **Configuration** dropdown is always active, even after a sample from the Experiment has been recorded.

## Update all samples

**Update all samples** applies the change of target and label names to all Samples in an Experiment or Group (instead of only to the current sample, which is the default).



- Click **Update all samples**, then select from the following options in the dropdown list:
  - **Apply to Experiment**
  - **Apply to Group**
- **Apply to Experiment:** Applies the changes to all Samples in the Experiment. This option keeps the IS/badge state of all existing samples unchanged as the Experiment-level IS are updated.
- **Apply to Group:** Applies the changes to all Samples in the Group and updates the IS/badge state of all existing Samples according to the following criteria:
  - Any Sample with an IS badge remains unchanged (i.e., the badge remains)
  - If the Auto Update Experiment IS is checked and the active Sample does not have data, any Sample in the Group that does not have an IS badge remains unchanged (i.e., no IS badge).
  - If the Auto Update Experiment IS is unchecked or the active Sample has data, any Sample that did not have an IS badge is assigned an IS badge.
- The **Update all samples** option is inactive during acquisition and when processing a Plate. When acquisition is paused, the button is active.

## Parameter selector

**Parameter selector** checkboxes allow you to select parameters (including **Event** number) and their respective measurements for any channel available in the selected configuration.

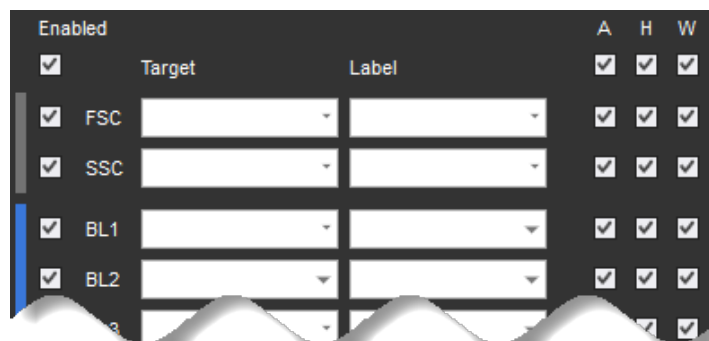
Enabled	Target	Label	A	H	W
<input checked="" type="checkbox"/>	FSC		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	SSC		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	BL1		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	BL2		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	BL3		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	RL1		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	RL2		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	RL3		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	VL1		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	VL2		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	VL3		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	VL4		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	YL1		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	YL2		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	YL3		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
<input checked="" type="checkbox"/>	YL4		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>

- By default, all parameters are selected except for the **Event** number parameter.
- Selecting the **Enable** checkbox enables all available channels with a single click.
- **Area**, **Height**, and **Width** measurements can be enabled for all channels with a single click by selecting the respective single-click checkbox (**A**, **H**, **W**).
- If all parameters or measurements are enabled using a single-click checkbox, then all related checkboxes are selected in that column.  
If any of those checkboxes are deselected, then the respective single-click checkbox is unchecked.

- Deselecting a **channel selector** checkbox unchecks all measurement checkboxes for that channel and disables the **Target** and **Label** fields.  
The **channel selector** checkbox is selected when at least one measurement is checked for that channel.  
If all measurements for a channel are unchecked, the **channel selector** checkbox is unselected and the **Target** and **Label** fields are inactive.
- **Time** parameter cannot be deselected.
- Parameters cannot be enabled or disabled in the current sample (wells or tubes) after that sample has been recorded and contains an FCS file.
- Parameters can be enabled or disabled during acquisition and before recording a sample. When parameters are enabled or disabled, the data will be cleared.
- Any parameter that is used as a Compensation control is disabled.

## Target textbox

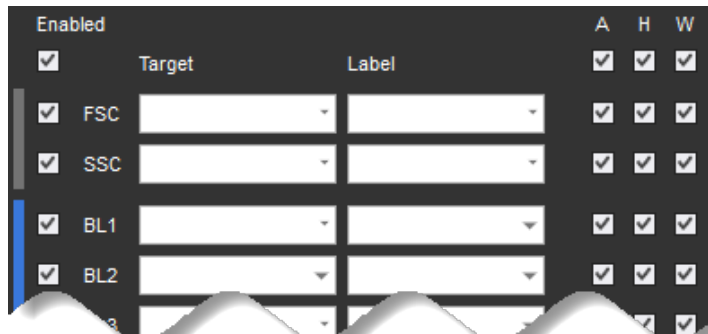
- **Target** textbox enables you to enter the target name of the label in the text field or leave it blank. The **Target** field is available for scatter and fluorescence parameters only.



- You can update the target name before or after recording.
- Any changes made in the **Target** field updates the parameter name used in the Workspace.
- Entries made in the Target textbox are used by the **Smart Gate Naming** tool as described in “Name options” on page 459.
- You can enter up to 50 characters into the **Target** textbox. Characters above this limit cannot be entered. Any ASCII character is valid.  
The tooltip for Target textbox reads “Example: CD3”.
- Any changes made to the target name post-recording only applies to the active sample, unless **Update all samples** button is clicked, which applies the target name to all samples in the Experiment.
- The tooltip for the **Update all samples** button displays “Updates Target and Label names for all samples within experiment.”

## Label textbox

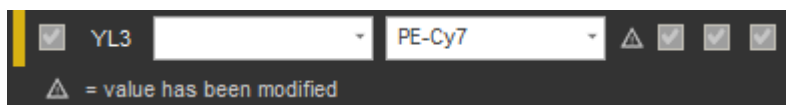
- The **Label** textbox enables you to enter the fluorophore name (i.e., label name) in the text field or leave it blank. The **Label** field is available for scatter and fluorescence parameters only.



- You can update the **Label** name before or after recording.
- You can enter up to 50 characters into the textbox. Characters above this limit cannot be entered. Any ASCII character is valid. The tooltip for **Label** textbox reads “Example: FITC”.
- Any changes made to the **Label** name after recording only applies to the current sample, unless **Update all samples** button is pressed, which applies the fluorophore name to all Samples in the Experiment.

## Notification badge

If you change the **Configuration**, **Target**, or **Label** for any sample after recording, a **notification badge** is displayed next to adjacent to the affected parameter or the instrument **Configuration** dropdown (if changed).

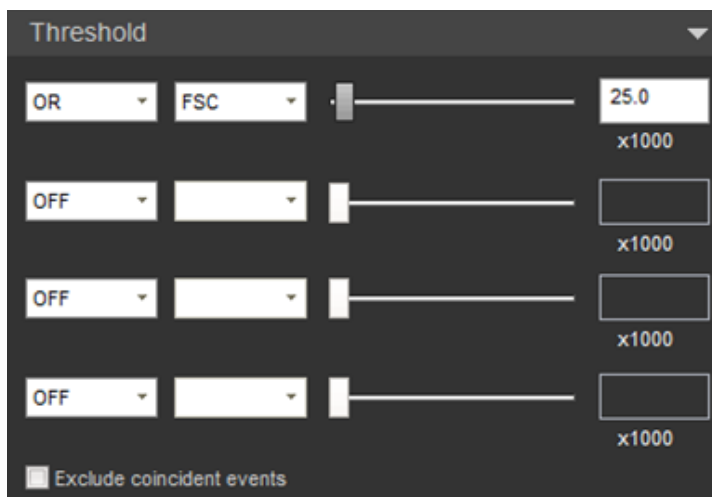


- Clicking the **notification badge** restores the **Target** or **Label** to the original value stored in the FCS file.
- Clicking the button next to the **Configuration** combo box restores the selection to the configuration used at the time the file was recorded.

# Threshold

## Overview

*Threshold* settings allow you to set the minimum signal level for up to four channels (i.e., detector) to eliminate unwanted events and reduce noise. You can also combine the thresholds using Boolean operators (see “Logic selector” on page 393).



- Threshold settings contain the following controls in the order listed:
  - *Logic selector combo box* (“Logic selector” on page 393)
  - *Channel selector combo box* (“Channel selector” on page 393)
  - *Threshold adjustment slider* (“Threshold adjustment slider bar” on page 394)
  - *Threshold signal textbox* (“Threshold textbox” on page 394)
  - *Exclude coincident events checkbox* (“Exclude coincident events check box” on page 395)
- A new value for the threshold signal is applied when the slider bar is moved, if the threshold logic or channel is changed, when the focus is lost from the threshold textbox, or when the **Enter** key is pressed.
- When a new threshold value is applied, existing data are cleared, and the software records and analyzes only the events with parameter values above the set threshold.
- Adjustments made to threshold with the slider bars dynamically refresh the data on the Workspace.

## Logic selector

**Logic selector** allows you to select up to four channels to apply threshold(s) and to combine the thresholds using Boolean operators.



- Each logic selector dropdown includes three options: **AND**, **OR**, and **OFF**.
- If the **OFF** option is selected for a channel, the detector is ignored for threshold purposes, and the channel selector, slider, and textbox are grayed out.
- The **AND** and **OR** selections follow Boolean logic rules.
  - If the **AND** operator is selected, all the set thresholds must be met before the data is collected.
  - If the **OR** operator is selected, data collection begins when at least one of the thresholds is met.
- Since the selected logic operator determines the enabled/disabled state for the rest of the controls, the Logic selector combo boxes are available before the other threshold controls.

## Channel selector

**Channel selector** allows you to select the channels to apply threshold.



- The first dropdown list includes all enabled channels selected in the Parameters group (excluding Time and Event) (“Parameter selector” on page 389).
- Each subsequent dropdown list shows only the remaining available parameters. For example, if FSC is used as the first threshold, then all other dropdowns will not have FSC as an option.
- Threshold is only applied to parameters that are visible, allowing up to four threshold permutations.
- Default Threshold settings are: **OR** (selected in the first Logic combo box), **FSC** with a value of 25 (× 1000), and **OFF** on the second, third, and fourth logic dropdowns (i.e., not used for threshold).
- When other Threshold settings are added, their default value will be 10 (× 1000).

## Threshold adjustment slider bar

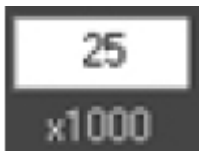
**Threshold adjustment slider bar** allows you to adjust the numerical threshold setting.



- Adjusting this slider dynamically updates the value in the threshold textbox accordingly. While adjusting the slider bar, a tooltip displays the full threshold value (i.e., the textbox value  $\times$  1000).
- The lower limit is 0.1. The upper limit is 1/1000 of the range set for the height measurement of the selected channel.
- Moving the slider bar adjusts the threshold setting by increments of 10. When starting at the lower limit of 0.1, the next level will be 10, after which the threshold will increase in steps of 10.
- You can use the **up/right** and **down/left arrow** keys can also be used to adjust the slider bar in 1 unit ( $\times$  1000) increments. The up/right arrow keys increase the value, and the down/left arrow keys decrease the value.
- When the textbox or slider bar is in focus, you can use the **mouse scroll wheel** to adjust the slider bar position.

## Threshold textbox

**Threshold** textbox allows you to type in the numeric value for a threshold setting. The software automatically multiplies the input value by 1000.



- Only numeric values are allowed in the textbox. One decimal place is always displayed. If you do not enter a decimal place, it will be automatically added with a zero in the tenths place.
- Changing the value in the textbox automatically updates the position of the slider after the textbox loses focus, the **Enter** key is pressed, the **up/right** and **down/left arrow** keys are pressed, or when the value is changed using the **mouse scroll wheel**.
- The accepted range of values is 0.1 for the lower limit, and 1/1000 of the allowed maximum for the height measurement for the upper limit (220 by default). The resolution of entry is 0.1.
- When a number is entered that is greater than the maximum allowed value for the parameter, the focus is set to the entire text to allow editing (instead of defaulting to maximum).
- If the values are deleted from the textbox and the textbox loses focus, then the previous value is restored to the textbox.



## Exclude coincident events check box

**Exclude coincident events** check box removes coincident events.

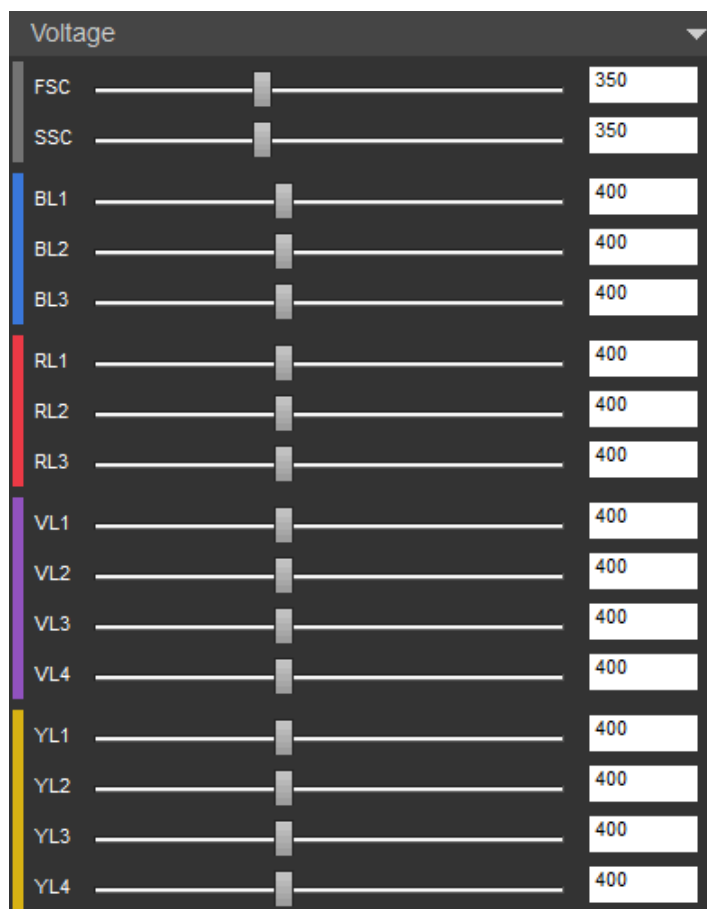


- When checked, the coincident events are removed from the data stream, and the number of aborted coincident events is counted in the FCS file. Aborted coincident events are annotated using the \$ABRT keyword.
- When checked, the number of aborted coincident events is updated in real-time in the Collection Panel.
- By default the checkbox is unchecked.

## Voltage

### Overview

**Voltage** settings allow you to adjust voltages for each channel. The number of channels shown depends on the configuration of the instrument.



- **Voltage** settings has the following controls:
  - **Voltage adjustment slider** (“Voltage adjustment slider bar” on page 397)
  - **Voltage value** textbox (“Voltage value textbox” on page 397)
- By default, FSC and SSC voltages are set at 350 and the fluorescent parameters are set at 400. Upon accepting the recommendations of BFR (“Baseline Functional Response (BFR)” on page 540), the voltages are updated with the optimal PMTVs.
- You can adjust the PMT voltage for an available channel using the Voltage adjustment slider or the textbox for that channel.  
If a parameter is deselected in the Parameters group (“Parameter selector” on page 389), then the deselected channel is grayed out.
- Adjustments made to voltages with the Voltage slider dynamically update the data on the Workspace. However, adjustments are not applied until the user stops pressing the mouse button when using the Voltage slider.

- Adjustments made using the textboxes automatically update the data on the Workspace when you click elsewhere on the screen or press the **Enter** key.
- When Compensation samples are run, the respective channel is highlighted with a light blue color.
- When a compensation channel is recorded or the Compensation sample contains an FCS file, all controls for other fluorescence channels are inactivated for the remaining samples in the Experiment.

---

**Note:** This ensures that the voltages used in the compensation calculation are not modified and the parameters are not inactivated.

---

- If the compensation settings are removed from the Experiment, the parameter and voltage controls for the fluorescence channels are reactivated.

## Voltage adjustment slider bar

**Voltage adjustment slider bar** allows you to adjust the PMT voltages for available channels.



- Adjusting the voltage slider bar dynamically updates the value in the voltage textbox accordingly. However, adjustments are not applied until the user stops pressing the mouse button when using the slider.
- The range of the voltage slider bar is from 1 to 1000 (left to right).
- Moving the voltage slider bar adjusts the PMT voltage by increments of 20.
- When the voltage slider or textbox is in focus, you can use the **mouse scroll wheel** or the **up/right** and **down/left arrow** keys to adjust the slider bar position by increments of 10. The up/right arrow keys increase the value, and the down/left arrow keys decrease the value.

## Voltage value textbox

**Voltage value** textbox allows you to directly enter the desired voltage setting for a specific channel.



- Only numeric values are allowed in the textbox. Non-numeric characters are not registered.
- The accepted range of values is 1–1000 mV for all channels. The resolution of entry is 1 mV.
- Changing the value in the textbox automatically updates the position of the slider after the textbox loses focus, the **Enter** key is pressed, the **up/right** and **down/left arrow** keys are pressed, or when the value is changed using the mouse scroll wheel.
- Adjustments to the textbox refresh the data on the Workspace once focus is lost or **Enter** is pressed.
- When a number is entered that is greater than the maximum allowed value for the parameter, the focus is set to the entire text to allow editing (instead of defaulting to maximum).

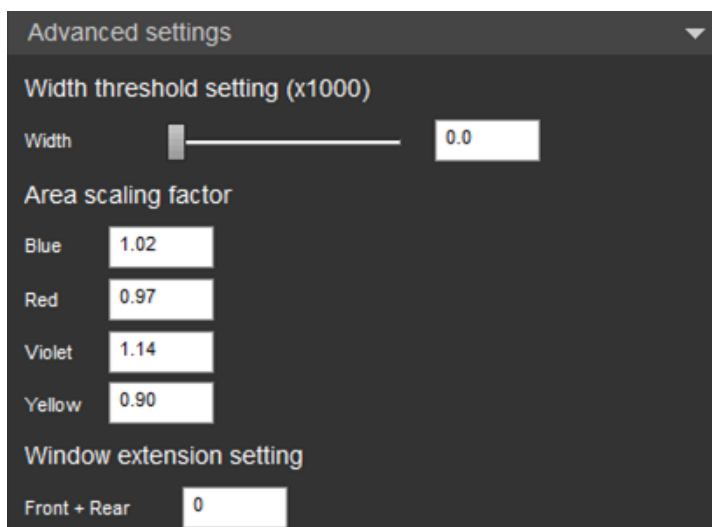
- If the values are deleted from the textbox and the textbox loses focus, then the previous value is restored to the textbox.
- During compensation, the fluorescence channels are highlighted with a light blue color.

## Advanced™ settings

### Overview

Advanced™ Instrument Settings allows you to adjust the following settings, if you have the appropriate permissions (see Note below):

- *Width threshold setting* (“Width threshold setting” on page 399)
- *Area scaling factors (ASF)* (“Area scaling factor (ASF)” on page 400)
- *Window extension setting* (“Window extension setting” on page 401)



- If you do not have permissions to any of the functions within this section, the Advanced™ settings section is not visible.
- Adjustments to the slider bar dynamically refresh the data on the Workspace. Adjustments to the textboxes update the data once you click elsewhere on the screen or when you press the **Enter** key.
- If the values are reset to default, the data is cleared and new values take effect.
- If you modify the area scaling factor (ASF) for any laser or change the window extension setting, a *notification button* (“Notification button” on page 401) is displayed next to the affected setting, and a notification key is shown at the bottom of the panel.
- Changes made to Advanced™ settings persist for the current user and the current experiment. New experiments will use the default system settings.
- Changes made to Advanced™ settings for one user are not applied to other users.

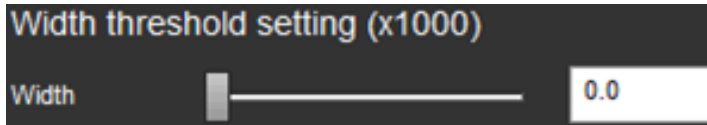
---

**Note:** The Advanced™ settings options are available only to users with Advanced™ User and Administrator level permissions. For more information, see “Account permissions” on page 44.

---

## Width threshold setting

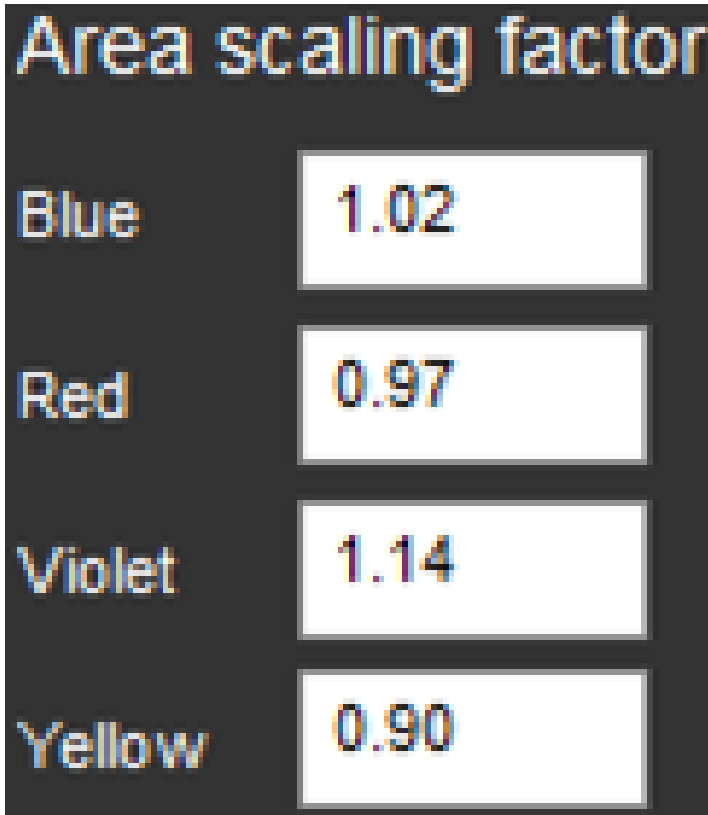
**Width threshold setting** sets the threshold above which the width measurement is determined. It consists of a slider and a textbox.



- The width threshold setting is set per Experiment. When it is saved as default instrument settings, it is used for new Experiments.
- The range of the width threshold setting is from 01 to 1048.5 (left to right), with a resolution of 0.1. The entered values are in thousands. The default value is 1.
- Moving the slider bar adjusts the value by increments of 10.
- Only numeric values are allowed in the textbox. One decimal place is always displayed. If you do not enter a decimal place, it is automatically added with a zero in the tenths place.
- When the slider or text box is in focus, you can use the **mouse slider wheel** or the **up/right** and **down/left arrow** keys to adjust the slider bar position by increments of 5.
- Changing the value in the textbox automatically updates the position of the slider after the textbox loses focus, the **Enter** key is pressed, the **up/right** and **down/left arrow** keys are pressed, or when the value is changed using the mouse scroll wheel.
- If the values are deleted from the textbox and the textbox loses focus, then the previous value is restored to the textbox.

## Area scaling factor (ASF)

**Area scaling factor (ASF)** option sets the value to scale height and area to equivalent values. Area scaling is calculated during the Performance Test and is automatically applied to new Experiments created after Performance Test is completed.

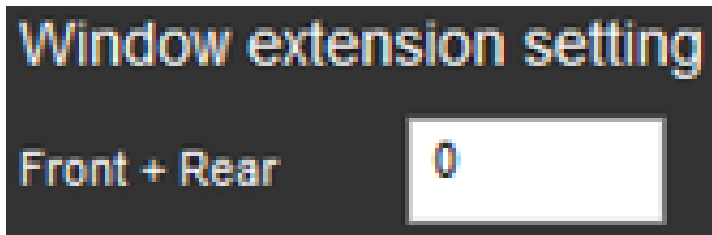


Area scaling factor	
Blue	1.02
Red	0.97
Violet	1.14
Yellow	0.90

- The ASF can be set per Experiment or it can be used for all new Experiments when saved as default Instrument Settings.
- The ASF can be set for each configured laser and only configured lasers are listed.
- The ASF values are updated after each successful Performance Test. The default values are set by the most recent, successful Performance Test. If no Performance Test or Baseline has been run, the default values are set to 1.0
- The ASF values range from 0.1 to 10.
- Only numeric values are allowed in the textbox. Non-numeric characters are not registered.
- If the System Settings update the Area scaling factor settings, the ASF values are updated to the System Settings.
- If the ASF value is modified in the text box for a specific laser, the System Setting is overridden such that the ASF value for that modified laser will not automatically update based on each successful Performance Test.
- If the ASF values are deleted from the textbox and the textbox loses focus, the previous value is restored to the textbox.
- If you do not have permission to access the Advanced™ Instrument Settings, the ASF value used is set by the System Instrument Settings.

## Window extension setting

The **Front + Rear window extensions** can be configured for an Experiment and saved as part of the default Instrument Settings.



- The value ranges from –40 to 75. The default value is 0.
- Only integer values are allowed in this text box.
- The window extension value input in the text box corresponds to a single extension and the full added extension is 2X of the input value.
- If the value is deleted from the text box, and the text box loses focus then the previous value will be restored to the text box.
- If the value is modified, the system window extension setting is overridden such that the value will not automatically update based on each successful Performance Test.

## Notification button

If you modify the area scaling factor (ASF) value for any laser or change the window extension setting, a *notification badge* is displayed adjacent to the affected setting.

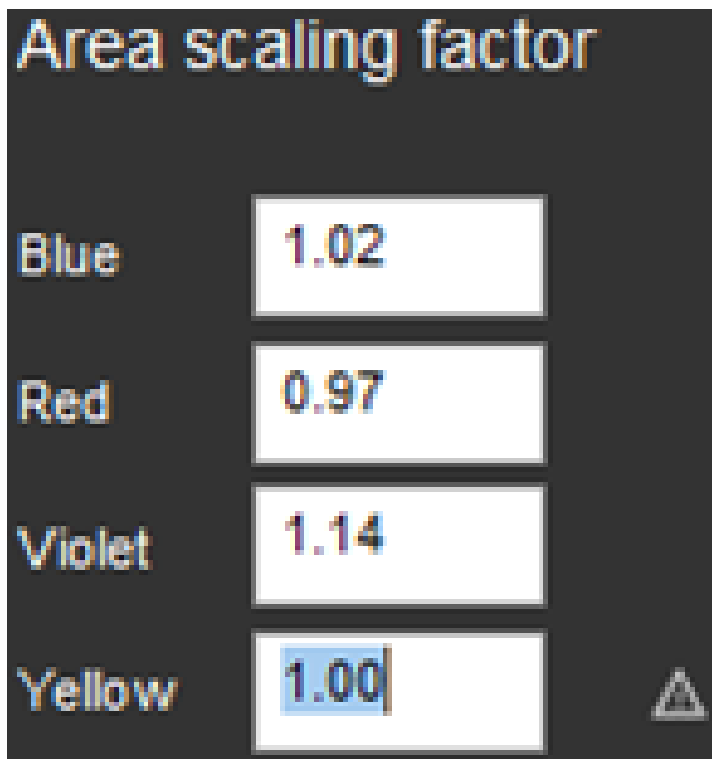
- Clicking the notification button next to the area scaling factor (ASF) restores the affected setting to the System Setting value for the corresponding laser.
- Clicking the notification button next to the window extension setting restores the setting to the System Setting value.

## System instrument settings

### Overview

*System Instrument Settings* allows you to adjust the following system settings, if you have the appropriate permissions (see Note below):

- *Area scaling factor (ASF)* (“Area scaling factor (ASF)” on page 400)
- *Window extension setting* (“Window extension setting” on page 401)



- If you do not have permissions to any of the functions within this section, the System Instrument Settings section is not visible.
- If the values are reset to default, the data is cleared and new values take effect.
- If you modify the Area scaling factor (ASF) value for any laser or change the window extension setting, a *notification badge* is displayed adjacent to the affected setting, and a notification key becomes visible at the bottom of the panel (“Area scaling factor (ASF)” on page 400).
- Changes made to System Settings are applied to each new Experiment created after the values have been changed. These changes persist until the next Performance Test is completed or until the System Settings Values are changed by an authorized user.
- Changes made to System Settings are applied as default to all user accounts after the values are changed.
- The System settings are updated after each successful Performance Test.

---

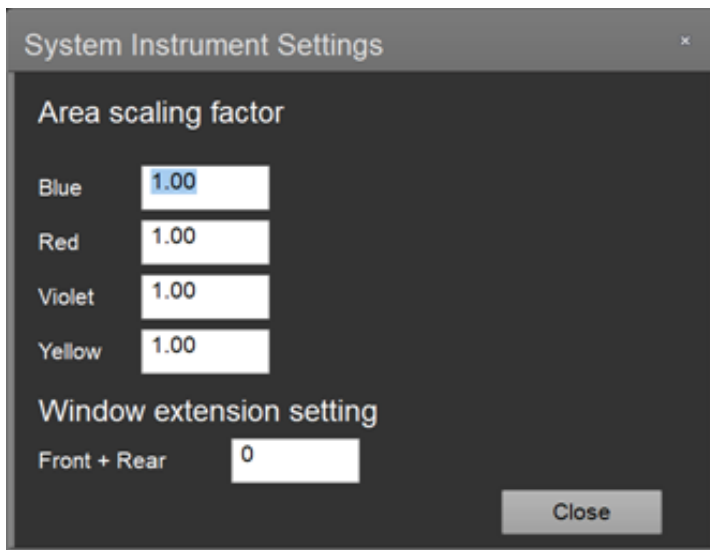
**Note:** The System Instrument Settings options are available only to users with Administrator level permissions. For more information, see “Account permissions” on page 44.

---



## Area scaling factor (ASF)

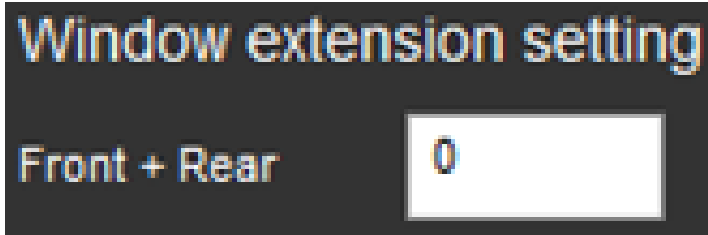
**Area scaling factor (ASF)** option sets the value to scale height and area to equivalent values. Area scaling is calculated during the Performance Test.



- The ASF setting applies to all Experiments and Users, where the Advanced™ Setting ASF value has not been set (“Area scaling factor (ASF)” on page 400).
- The ASF can be set for each configured laser and only configured lasers are listed.
- The ASF values are updated after each successful Performance Test. The default values are set by the most recent, successful Performance Test.
- If no Performance Test or Baseline has been run, the default values are set to 1.
- The values range from 0.1 to 10.
- Only numeric values are allowed in the textbox. Non-numeric characters are not registered.
- If the values are deleted from the textbox and the textbox loses focus, then the previous value is restored to the textbox.

## Window extension setting

The **Front + Rear window extensions** can be configured for an Experiment and saved as part of the default instrument settings.



- The value ranges from –40 to 75. The default value is 0.
- Only integer values will be allowed in this text box.
- The window extension value input in the text box corresponds to a single extension and the full added extension is 2X of the input value.
- If the value is deleted from the text box, and the text box loses focus then the previous value will be restored to the text box.
- The window extension value is reset to 0 after each successful Performance Test.

## Notification icon

If you modify the Area scaling factor (ASF) value for any laser or change the Window extension setting, a *notification badge* is displayed adjacent to the affected setting.

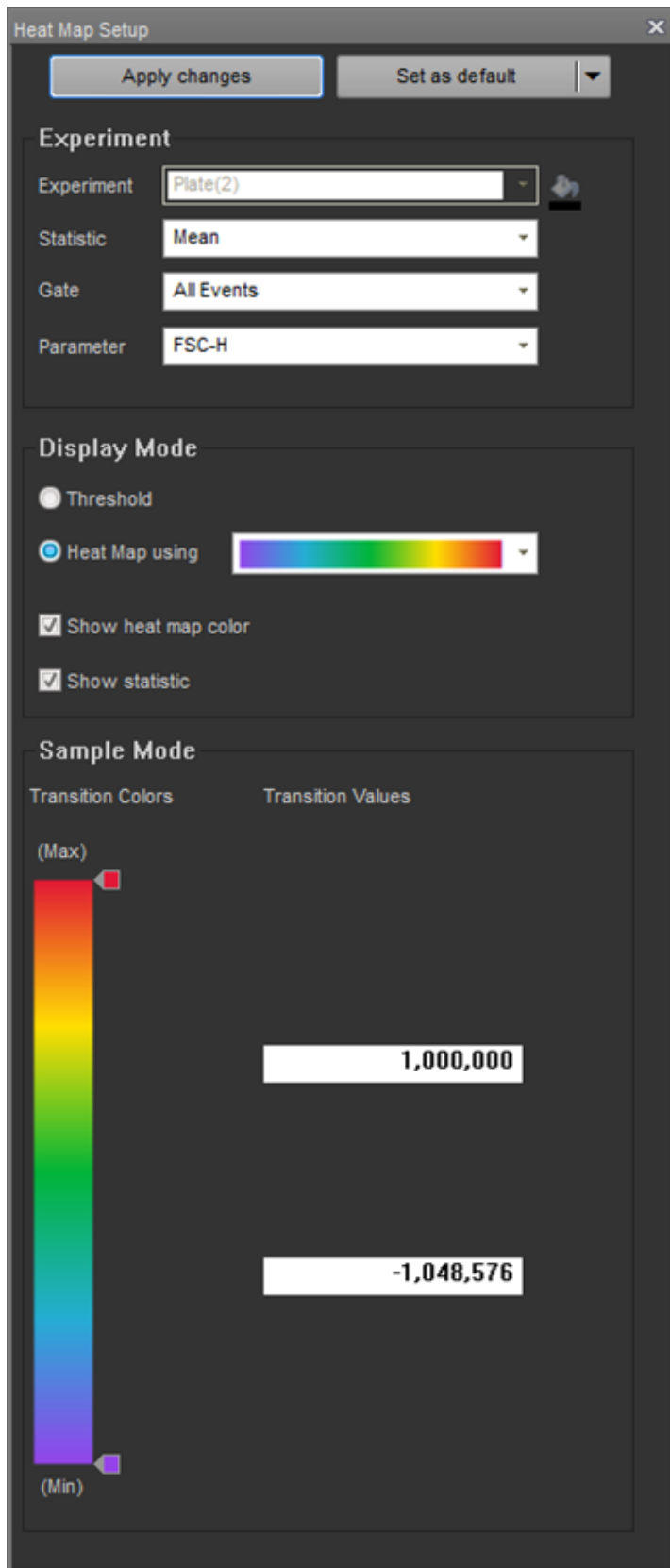
- Clicking the notification button next to the Area scaling factor (ASF) restores the affected setting to the most recent successful Performance Test value.
- Clicking the notification button next to the Window extension setting restores the setting to its default value of 0.



# Heat map setup panel

## Overview

*Heat Map Setup* panel allows you to select the statistic, gate, and parameter for visualizing the data from the current Experiment in the *Heat Map view* (Chapter 6, “Heat Map View”), and to specify the **display mode** and **transition values** to help you analyze the data at a glance.



- In the **Threshold display mode**, each well and/or tube is assigned a color based on whether the selected statistic for that sample falls above or below the set threshold value.
- In the **Heat Map display mode**, individual values in each well and/or tube are represented as colors on a gradient.
- By default, the Heat Map Setup panel is docked to the left of the Main Application Area. The docking properties of the panel are described on “Docking locations” on page 58.
- The Heat Map Setup panel is organized into four functional groups in the following order:
  - Apply, Save, and Load options (“Apply, save, and load options” on page 407)
  - Experiment (“Experiment” on page 408)
  - Display mode (“Display mode” on page 412)
  - Sample value (“Sample mode – threshold” on page 414)
- Sample value options on the Heat Map Setup panel are contextual; the controls displayed depend on whether you select Threshold or the Heat Map in the Display mode options (“Display mode options” on page 412).
- By default, the Heat Map Setup panel displays the sample value controls for the Threshold mode.

---

**Note:** For information on the Heat Map view, including Sample display and Experiment Setup, refer to “Heat Map View” on Chapter 6, “Heat Map View”.

---

## Apply, save, and load options

### Overview

*Apply, Save, and Load options* consist of the two controls that lie at the top of the Heat Map Setup panel.



### Apply changes

- **Apply changes** applies the current Heat Map settings to all samples in the current Experiment.
- This option is active if the Heat Map settings of the current Experiment have been changed.
- If the Heat Map Setup panel loses focus after a change has been made, the new settings are automatically applied to the Heat Map.

The Status Notification Display located above the Main Application Workspace (“Main application workspace” on page 56) then displays “The Heat Map settings have been updated” and provides an **Undo** option.

## Set as default split button

**Set as default** button is a split button that allows you to apply the Heat Map settings to the current Experiment, set them as default, or save (export) and retrieve them for other Experiments.

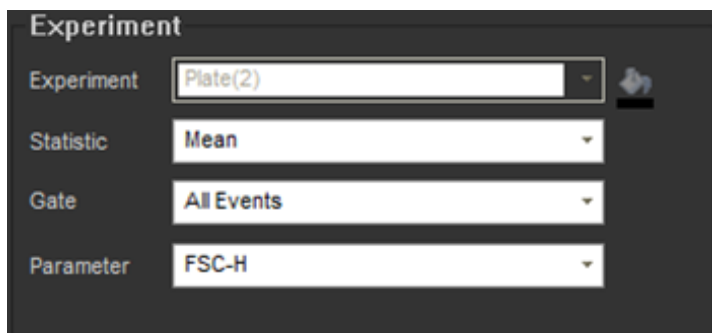
The main button on the left is the **Set as default** command and the right side is a dropdown containing **Load** and **Export** options.

- **Set as default:** Saves the current Heat Map setting as the default Heat Map settings for all new Experiments.
- **Load:** Opens the *File Browser dialog* (“Folder Browser dialog” on page 723).  
Using the File Browser dialog, you can select a Heat Map settings file from within any experiment in your folder, from available Shared folders, or from available default instrument settings files.
- **Export:** Opens the *File Save (Export) dialog* (“File Save (Export) dialog” on page 715).  
Using the File Save (Export) dialog, you can save the current Heat Map settings with a user-defined name in your own folder or any Shared folders in the database.  
You will be prompted for confirmation before overwriting an existing Heat Map settings file.

## Experiment

### Overview

*Experiment options* allow you to select the statistics, populations (all events or specific gated populations), and the parameter for the current Experiment to display in the Heat Map view. It also contains the control for assigning a unique indicator color for the current Experiment.



## Experiment dropdown

**Experiment** dropdown list displays the name of the Experiment defined within the current plate and any associated tubes.



- The Experiment dropdown control is disabled as each plate is limited to a single Experiment.
- The control is labeled with the name of the currently open Experiment. The default label is **Experiment**.
- Default settings are loaded the first time the Experiment Heat Map is selected.

## Experiment color button

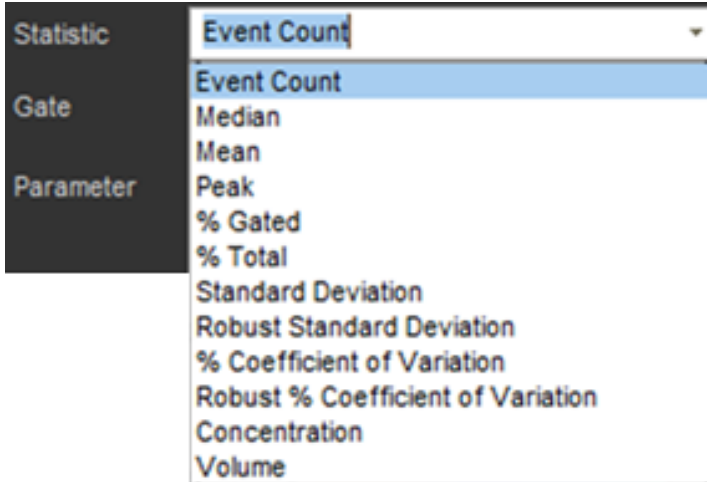
**Experiment color** button opens the *color picker dialog*, which allows you to select the indicator color for the current Experiment.



- By default, the Experiment color dropdown button shows the currently selected color for the Experiment.
- When applied, the selected color is reflected in the Heat Map view (Chapter 6, “Heat Map View”).

## Statistic dropdown

**Statistic** dropdown displays the list of available statistics and allows you to select the statistic of interest to display in the Heat Map view.

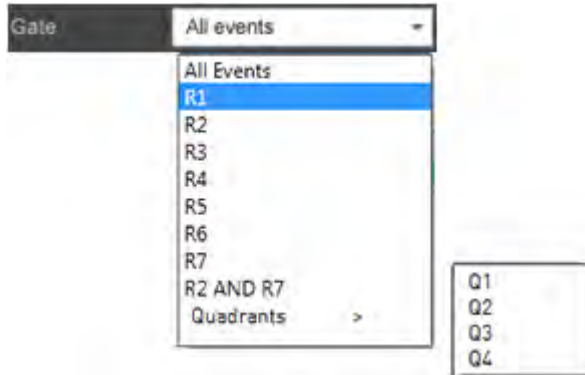


- The statistic dropdown control is labeled with the name of the currently selected statistic. The default statistic is **Event Count**.
- The dropdown list includes the following statistics:
  - Event Count
  - Median
  - Mean
  - Peak
  - % Gated
  - % Total
  - Standard Deviation
  - Robust Standard Deviation
  - % Coefficient of Variation
  - Robust % Coefficient of Variation
  - Concentration
  - Volume



## Gate dropdown

**Gate** dropdown list allows you to select the populations of interest (all events or specific gated populations).



- The dropdown list contains all gates in the Experiment-level Workspace (“Workspace levels” on page 118), including quadrant and derived gates.
- The default population is **All Events**.
- Only gates from the Experiment-level Workspace are displayed.
- Selecting a gate sets the parent gate for the Heat Map display.
- If the default Heat Map or loaded Heat Map refer to a gate that is not available on the Experiment workspace, the gate selection defaults to All Events.

## Parameter dropdown

**Parameter** dropdown allows you to select the parameter to display in the Heat Map view for the current Experiment.



- The Parameter dropdown is available only for intensity related statistics, Mean, Median, Standard Deviation (SD), Robust SD, Coefficient of Variation (CV), and Robust CV. It is hidden when Event Count, Percentage of Parent, Percentage of Total, or Concentration is selected in the Statistic dropdown list (“Statistic dropdown” on page 410).
- The dropdown list contains only the parameters on the plot for the gate selected in the Gate dropdown.
- If **All Events** is selected in the Gate dropdown, the parameters are not filtered and all enabled parameters are displayed by default.
- If the gate is derived from a Histogram plot (i.e., a single parameter plot), then the parameter list only contains the parameter that is on the selected plot. Selecting the **More** displays all available parameters.
- For a dual-parameter plot, the parameter list is filtered to show the two parameters that are on the plot from which the gate is derived. Selecting the **More** option shows all parameters.

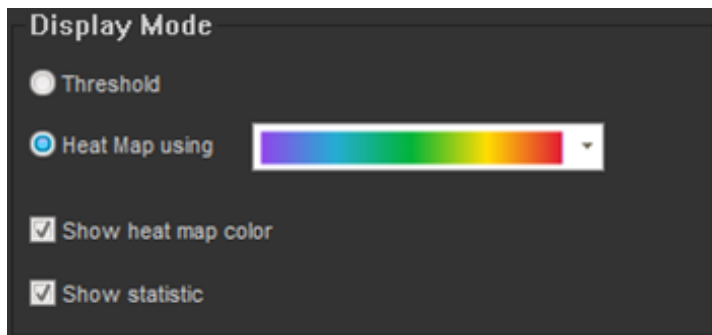
## Display mode

### Overview

*Display mode* options allow you to choose between the **Threshold** and **Heat Map** display modes and control the display of the colors used in the Heat Map view.

You can select the Threshold or the Heat Map mode by clicking the appropriate radio button.

The default option is Threshold mode.



---

**Note:** Threshold display mode is a visual aid to determine at a glance whether the signal from a well falls above or below a set value; it does not affect data collection. As such, it is different from the *Threshold settings*, which sets the minimum signal level for each detector to eliminate unwanted events and reduce noise. The threshold value set in Threshold settings instructs the software to record and analyze only the events with parameter values above the set threshold.

---

### Display mode options

#### Threshold mode

- **Threshold** radio button selects the Threshold mode and displays each well in user-defined colors based on whether the signal from the well and/or tube falls above or below the set threshold value.
- By default, only a single threshold is defined. However, you can create up to 10 signal threshold levels using the **Sample value** options (“Sample mode – threshold” on page 414).
- When **Threshold** mode is selected, choosing a color gradient from the dropdown switches the radio button to the **Heat Map** mode.

#### Heat map mode

- **Heat Map** radio button allows you to represent the data from each sample well as colors on a gradient selected from the **color gradient dropdown** list.

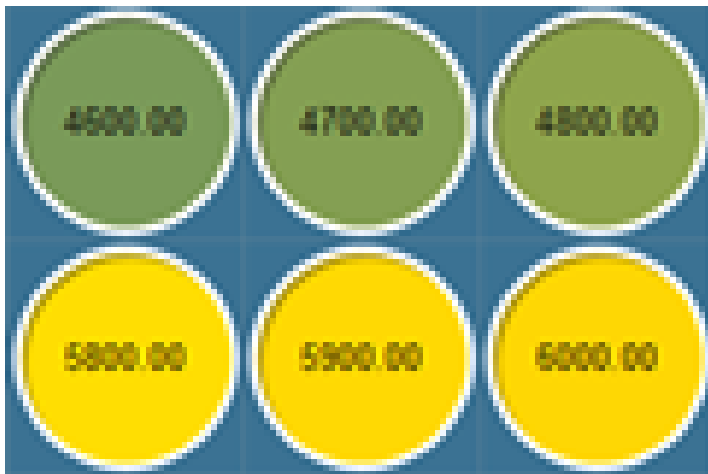


- There are six gradients available in the color gradient dropdown list, which are not editable. The default option is the first color range in the list.
- By default, only the minimum and maximum transition points are defined in Heat Map. However, you can create up to 10 transition points using the **Sample value** options (“Sample mode – threshold” on page 414).

### Show heat map color

- **Show heat map color** toggles the heat map color displayed on the Heat Map on and off.
- This option is selected by default (i.e., the Heat Map color is turned on).

### Show statistic

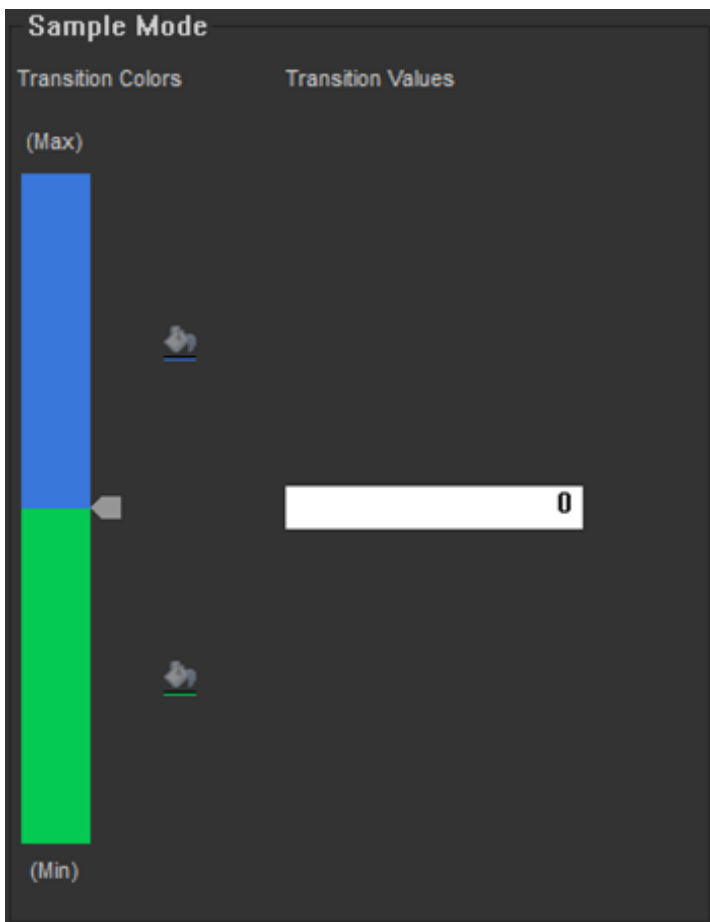


- When selected, the **Show statistic** option overlays the selected statistic centered in the wells and/or tubes.
- The background color includes the Experiment color (if selected) and the Heat Map color (if selected).

## Sample mode – threshold

### Overview

*Sample mode* options allow you to add and remove color transition points and set the data level and the associated color.



- Clicking the **Threshold** radio button in Display mode options (“Display mode options” on page 412) displays the sample value controls for the Threshold mode.
- In the Threshold mode, each well and/or tube is assigned a color based on whether the selected statistic for that sample falls above or below the set threshold value.
- Threshold mode is the default option for the sample value controls.

## Transition colors

- The color of the well in the Heat Map view (Chapter 6, “Heat Map View”) changes, if the threshold signal level is exceeded. It will not change if the value is the same or less than the transition value.
- By default, if the selected statistic is above the threshold, the well appears green; if it is below this level, the well appears blue.
- Clicking the **color picker** button next to a color block opens the standard color picker dialog, which allows you to define a different color for that block.



## Transition values

- By default, a single transition point (i.e., threshold) is set in the Threshold display mode. The default transition value is 10.
- You can enter the data value for a transition point directly in the **transition value** edit box.



- Transition value edit boxes accept only numerical characters. Decimal values are shown only if you enter them.
- Each transition value edit box allows manual entry of a number based on the statistic being displayed.
  - Percentage of count values: 0 to 100%.
  - Percent CV values:  $\pm 1000\%$
  - Events: 1 to 100,000,000
  - Intensity: minimum to maximum allowed for the selected parameter
- If the entry for the transition point exceeds the maximum value allowed, it will default to the maximum; if the value is lower than the minimum value allowed, then it will default to the minimum.
- If the statistic is changed, then the threshold returns to the default transition point set at 10.

## Create additional transition points

- You can add additional transition points by clicking on any given colored block sector. This inserts a new threshold bisecting the selected sector and adds an additional color block, which adopts the next color in the automatic color sequence.
- A transition value edit box is automatically created for each new level, and a color picker is associated with each new color block.
- You can create up to 10 levels. If you attempt to create more than 10 transition points, a warning message is displayed below the color bar.
- Clicking to create an additional transition point of a higher value creates a transition point at double the value of the first point.  
For example, if the first transition point is set at 100, the new one will be set to 200.
- Clicking to create an additional transition point of a lower value creates a transition point at half the value of the first point.  
For example, if the first transition point was set at 100, the new one will be set to 50.
- Clicking to create an additional transition point between two existing points creates a transition point half way between the two existing values.  
For example, if you click between the transition points set at 100 and 200, the new transition point value will be set to 150.
- Clicking and dragging an indicator arrow from the color bar by more than 20 pixels displays a red **X** icon, and when the mouse button is released, the selected transition point and the associated color picker and transition value edit box are removed.

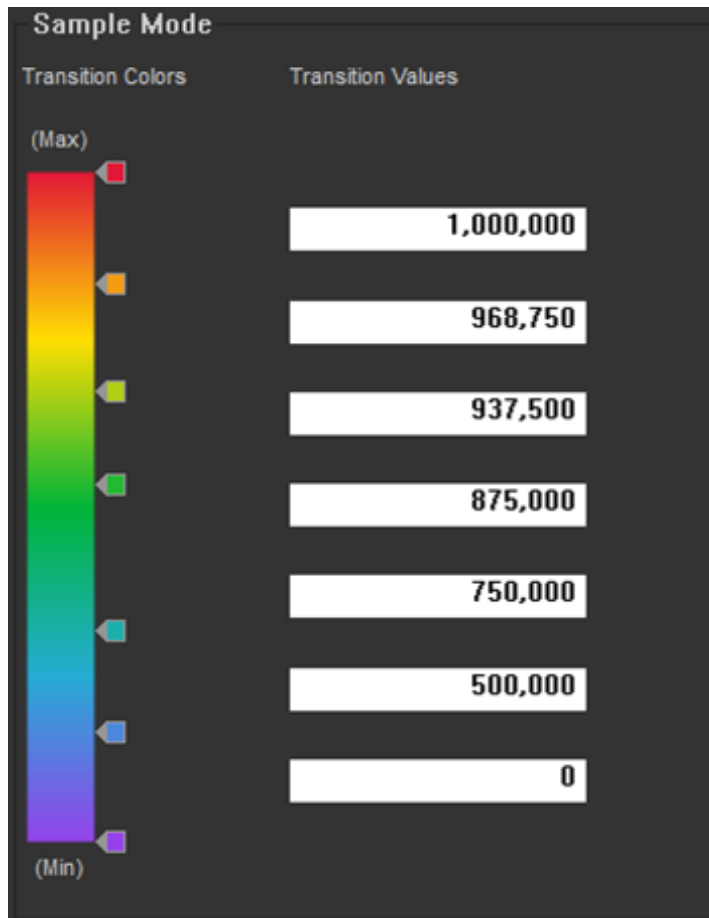


- There must always be one transition point defined. The last point cannot be removed.

## Sample mode – heat map

### Overview

*Sample mode* options allow you to add and remove color transition points and set the data level and the associated color.



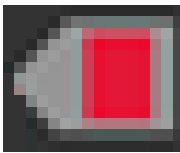
- Clicking the **Heat Map** radio button in Display mode options and selecting the transition colors from the **color gradient dropdown** list (“Display mode options” on page 412) displays the sample value controls for the Heat Map display mode (i.e., gradient mode).
- In the Heat Map mode, each well and/or tube displays the appropriate color from the selected gradient depending on whether the signal from the well falls within the set boundaries.

## Transition colors

- The color of the well in the Heat Map view (Chapter 6, “Heat Map View”) changes depending on where the result is in comparison to the selected color gradient.
- The Transition color gradient is selected using the **color gradient** dropdown list as described in “Display mode options” on page 412.



- Colored indicator arrows are displayed at the minimum and maximum points by default and any added transition points in between. The color displayed indicates the transition point color.



- Transition points can be moved up and down to adjust the color assigned to that transition point by clicking and dragging the indicator arrows with the mouse. The associated textbox will be highlighted in light blue to distinguish which transition point is being adjusted.

## Transition values

- By default, minimum and maximum transition points are defined in the Heat Map display mode.
- The default minimum transition point is 0 and the default maximum transition point is 10, 000, from bottom to top respectively. The default maximum for statistics containing a percentage is 100%.
- You can enter the data value for a transition point directly in the **transition value** edit box.
- Transition value edit boxes accept only numerical characters. Decimal values are shown only if you enter them.
- Each transition value edit box allows manual entry of a number based on the statistic being displayed.
  - Percentage of count values: 0 to 100%.
  - Percent CV values:  $\pm$  1000%
  - Events: 1 to 100,000,000
  - Intensity: minimum to maximum allowed for the selected parameter
- If the entry for the transition point exceeds the maximum value allowed, it will default to the maximum; if the value is lower than the minimum value allowed, then it will default to the minimum.
- If the statistic is changed, the minimum and maximum values revert to the appropriate default based on that statistic. Any middle transition points that existed are also removed.



## Create additional transition points

- You can add additional transition points by clicking on the color bar. This inserts a new transition point at the selected point and adds an additional indicator arrow, which adopts the color of the point selected.
- A transition value edit box is automatically created for each new transition point.
- You can create up to 10 levels. If you attempt to create more than 10 transition points, a warning message is displayed below the color bar.
- Clicking to create an additional transition point between two existing points creates a transition point half way between the two existing values.  
For example, if you click between the transition points set at 100 and 200, the new transition point value will be set to 150.
- The transition value entered must be in ascending order. If an incorrect value is entered, a warning balloon is displayed.
- Clicking and dragging an indicator arrow from the color bar by more than 20 pixels displays a red **X** icon, and when the mouse button is released, the selected transition point and the associated transition value edit box are removed.



- The minimum and maximum points cannot be removed.

## Overview

The *Customize panel* displays context sensitive formatting options relevant to objects selected on the Workspace or the Heat Map view.

Using the Customize panel, you can format the following objects:

- *Plots* (“Customize plot options” on page 421)
- *Text boxes* (“Customize text box options” on page 444)
- *Statistics boxes* (“Customize statistics box options” on page 445)
- *Gate styles and labeling* (“Customize gate options” on page 447)
- *Cell Image Container - Image Adjustment* (“Customize Image options” on page 484)

## Customize panel behavior

- To display the Customize panel, right-click an object to open its context menu, then select **Customize**.
- The formatting options in the Customize panel are context-sensitive and dynamically update depending on the currently selected objects.
- When a Compensation sample is active, all customize functionality is disabled except those specified in Chapter 21, “Compensation”.
- You can select multiple objects of the same type (e.g., plots), but not multiple objects of different types (e.g., plots and text boxes).

If you select multiple objects of different types, the Customize panel displays a warning message stating:

“Customize options are available only when similar objects are selected.”

# Customize plot options

## Overview

*Customize Plot options* allow you to customize properties for the following plot types: *Histogram*, *Dot*, *Density*, and *Precedence Density*.

---

**Note:** To set default options for plots, use the *Options dialog* (“Fonts and Styles” on page 649).

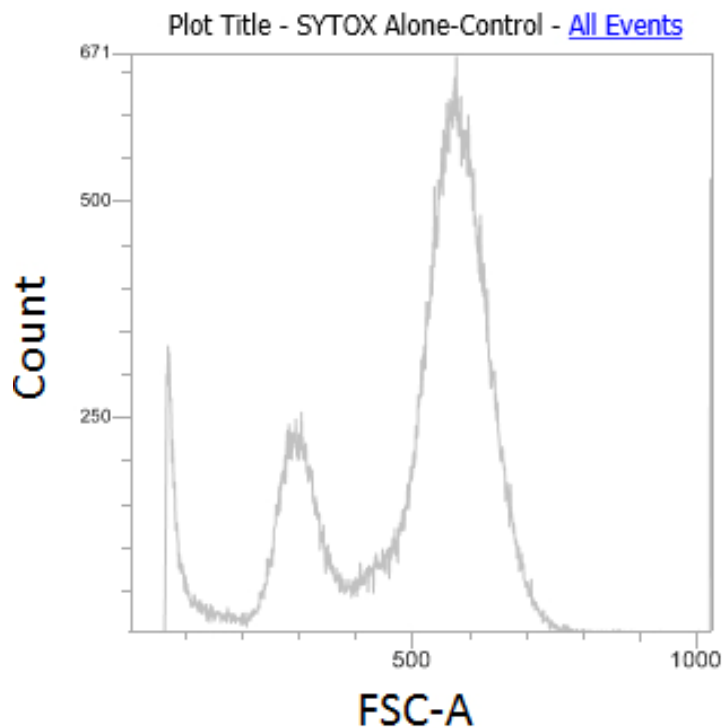
---

## Plot types

### Histogram plot

A *Histogram plot* is a graphical representation of single-parameter data and shows the relative number and distribution of events.

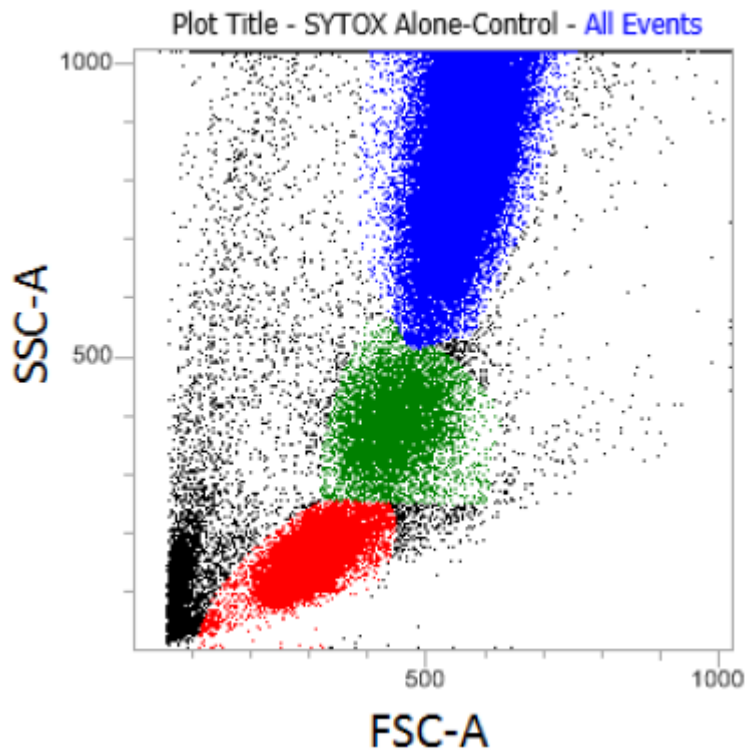
In a Histogram plot, the horizontal axis corresponds to the signal intensity of the selected parameter while the vertical axis represents the number of events (count).



## Dot plot

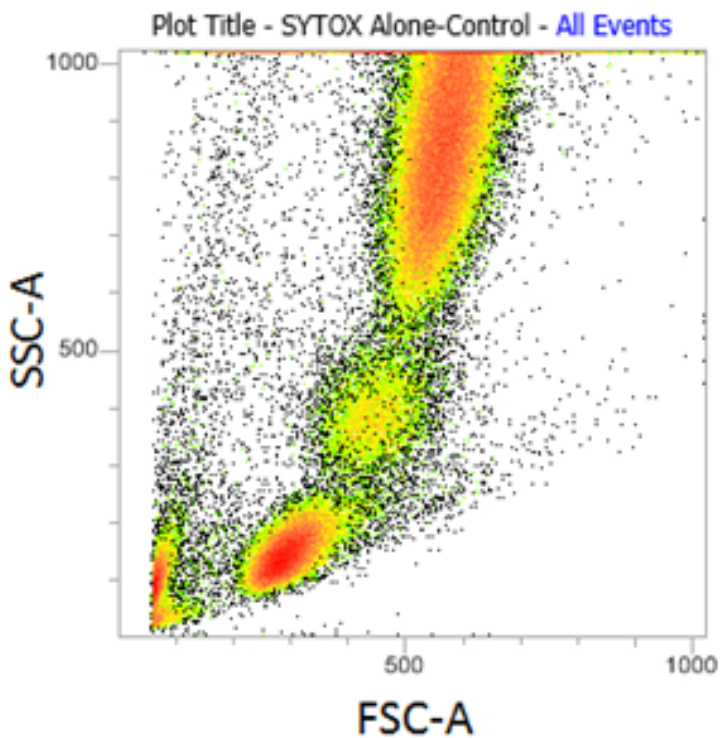
A *Dot plot* is a graphical representation of dual-parameter data, where each axis represents the signal intensity of one parameter.

Each dot in the plot corresponds to one or more events detected above the threshold. Different colors are used to represent the parent gate of events that fall within bins on the plots.



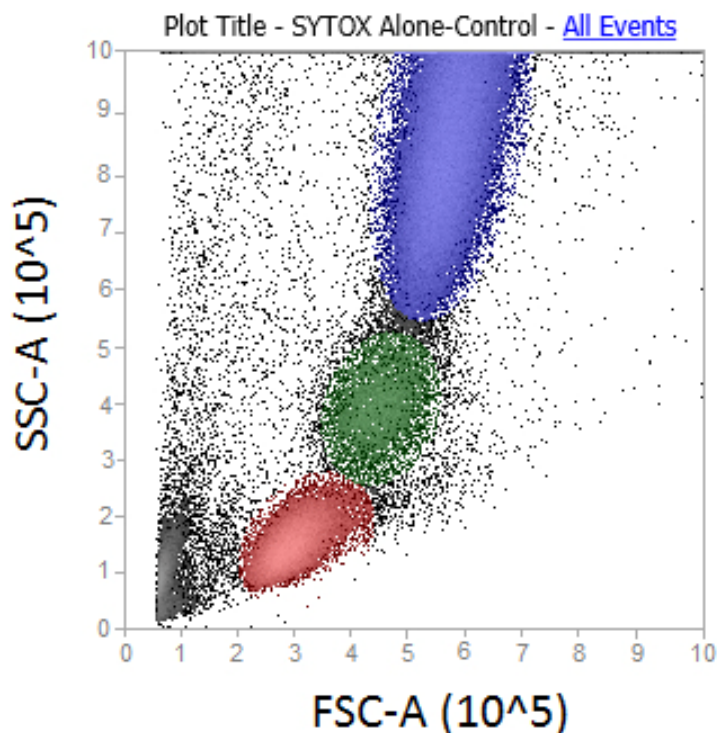
## Density plot

A *Density plot* is a graphical representation of dual-parameter data, where the colors represent the collection of events with the same intensity and each axis represents the signal intensity of one parameter.



## Precedence density plot

A *Precedence Density plot* is a combination of Dot and Density display, where a gradient is used to indicate the number of events within each of the plot bins and color is used to display the parent gate of events present.



## Customize panel behavior for plots

The Customize Plot options are displayed when one or more plots are selected. They are arranged into four groups:

- *General* (“General group” on page 427)
- *X-axis* (“X-axis group” on page 436)
- *Y-axis* (“Y-axis group” on page 441)
- *Text* (“Text group (plots only)” on page 442)
- The options displayed in each group vary, depending on the plot types and options selected. The example images below show the options available for Density plots when scatter or time and fluorescent parameters are selected.
- Each field in the panel displays the current settings for the selected plot. If multiple plots with differing values are selected, the value fields show the indeterminate state.

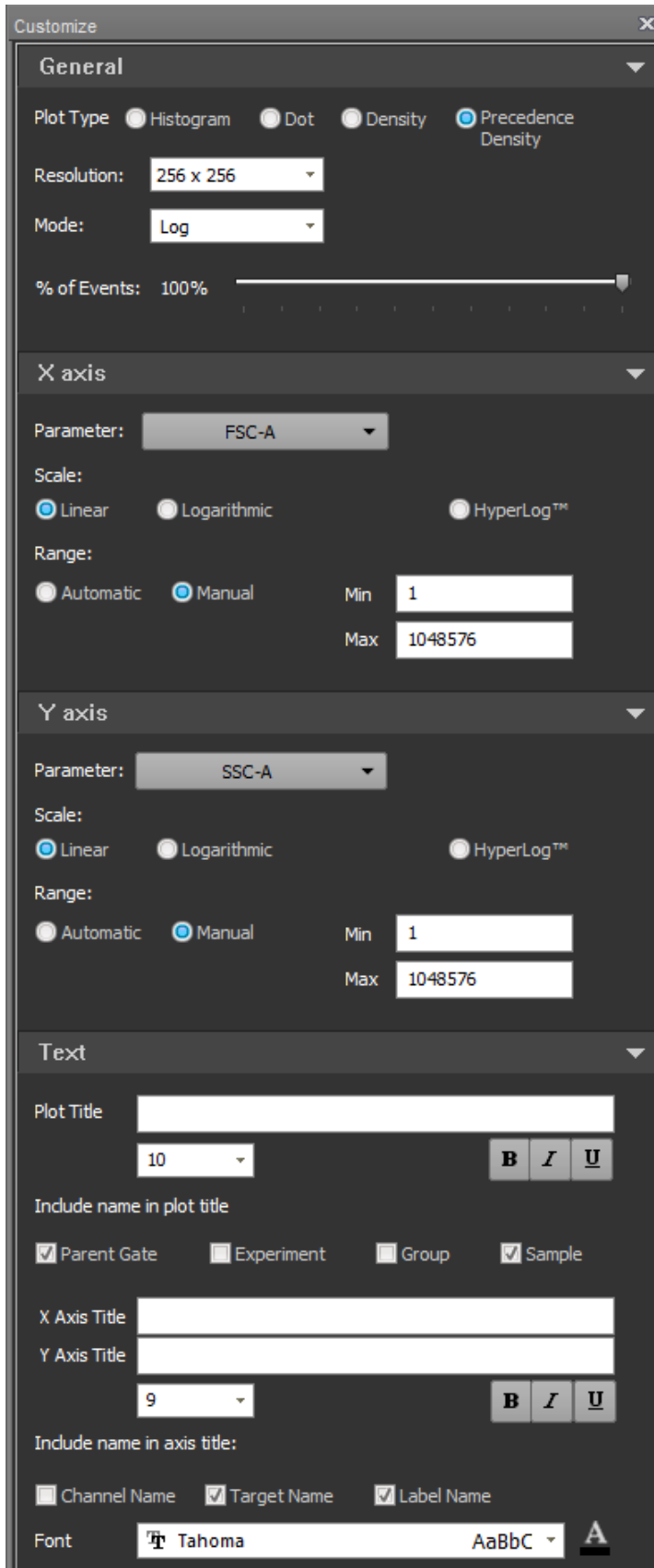
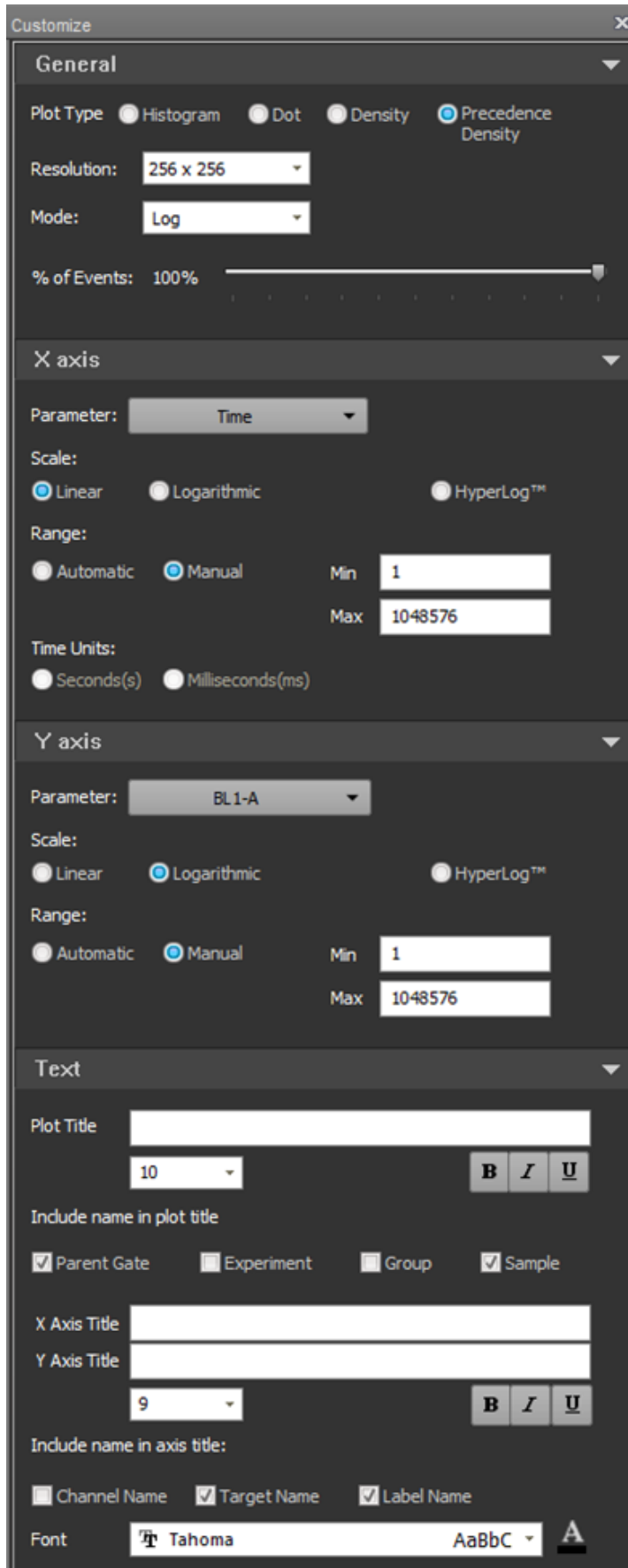


Figure 80 Customize Plot options for Density plots (scatter parameters)





## General group

*General group* controls the plot type and general plot characteristics (such as plot resolution).

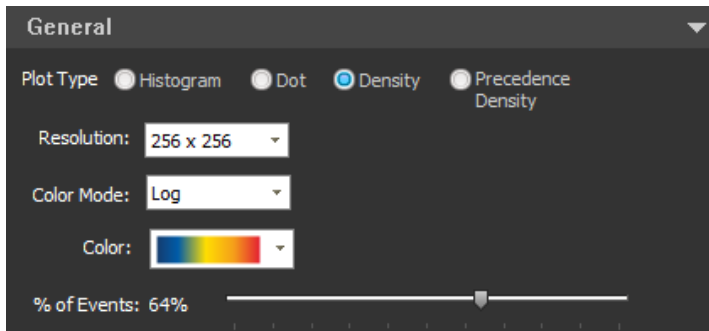
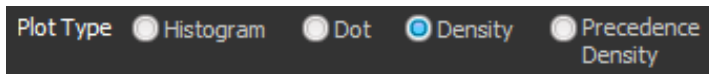


Figure 82 General group for Density plots

**Note:** The customization options that appear in the General group vary depending on the plot types selected. This section describes all possible options available in the General group. For illustrations of the plot-specific options, see “General group – plot-specific options” on page 434.

### Plot type radio buttons

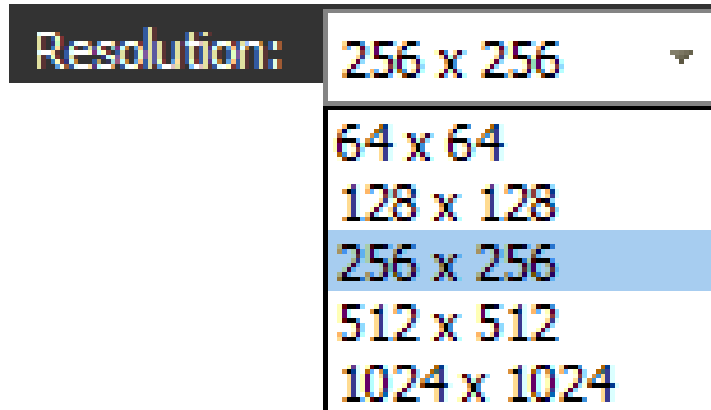
**Plot Type** radio buttons allow you to assign a plot type for the selected plots.



- The Plot Type option applies to all plot types and it is displayed when any plot is selected.
- If one or more plots of the same type are selected, the radio button for that plot type is selected.
- If multiple plot types are selected, all radio buttons are deselected.
- Selecting a radio button changes all selected plots to the selected plot type.
- If the current Workspace is a Compensation Workspace, the Histogram radio button is disabled when Density, Dot, or Precedence Density plots are selected. Similarly, if Histogram plots are selected, the Density, Dot, and Precedence Density radio buttons are disabled.

## Resolution dropdown menu

**Resolution** dropdown menu allows you to select a channel resolution for the selected plots.

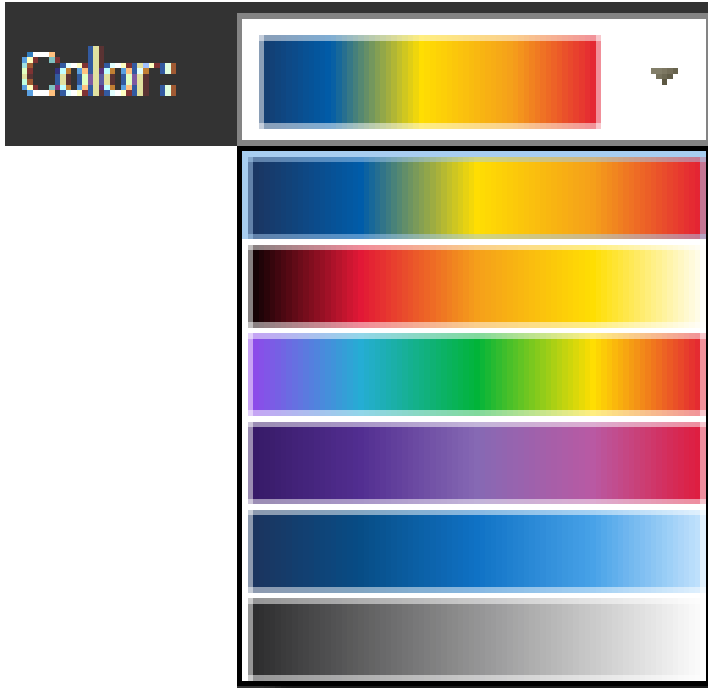


- Resolution option applies to all plot types and it is displayed when any plot is selected.
- If a single plot is selected, or if multiple plots with the same resolution (including a mixture of plot types) are selected, the Resolution dropdown menu displays the resolution of the currently selected plots.
- If multiple plots of any type with different resolutions are selected, the Resolution field is empty (blank).
- The maximum channel resolution is 1024 × 1024 channels for Dot, Density, and Precedence Density plots, and 1024 channels for Histogram plots.
- If a Histogram plot is selected along with two-parameter plots (Dot, Density, and/or Precedence Density plots), the Resolution dropdown menu for dual-parameter plots is displayed.
- If a Histogram plot is among the selected plots, the resolution is applied to the X-axis of the Histogram plot.
- The default resolutions are:
  - Histogram – 1024 channels
  - Dot – 256 × 256 channels
  - Density – 256 × 256 channels
  - Precedence Density – 256 × 256 channels

The defaults can be modified in *Plot Options* in the *Options dialog* (“Fonts and Styles” on page 649).

## Color dropdown menu

The **Color** dropdown menu allows you to select a color scheme for Density plots.



- Color option is displayed when one or more Density plots are selected.
- If a single Density plot is selected, or if multiple Density plots with the same color scheme are selected, the Color field displays the density color scheme of the currently selected plots.
- If multiple Density plots with different color schemes are selected, the Color field is empty (blank).
- You can select a different color scheme from the dropdown menu. The selected color is applied to all selected Density plots.
- The default color scheme can be modified in *Plot Options* in the *Options dialog* (“Fonts and Styles” on page 649).

## % of events slider

The **% of Events** slider adjusts the percentage of events displayed on the selected plots. The sample data is taken from the entire data file.



- % of Events option applies only to dual-parameter plots (Dot, Density, and Precedence Density plots), and is displayed when one or more dual-parameter plots are selected.
- % of Events option is not displayed when only Histogram plots are selected. If Histograms plots are selected along with two-parameter plots, adjusting the slider has no effect on the Histogram plots.
- The current value for the slider setting is shown to the left of the slider and it is updated as the slider is moved.
- When using the mouse, you can adjust in 1-unit increments.  
When using the arrow keys, you can adjust in 5-unit increments.  
Clicking the bar to the left or right of the slider moves the slider by 25 units.

## Color mode dropdown menu

**Color Mode** dropdown menu allows you choose **Linear** or **Log** binning of the data to the selected color palette for Density and Precedence Density plots.



- Color Mode option applies only to Density and Precedence Density plots and it is displayed when one or more Density and/or Precedence Density plots are selected.
- If a single Density or Precedence Density plot is selected, or if multiple Density and/or Precedence Density plots with the same color mode are selected, the Mode field displays the color mode of the currently selected plots.
- If multiple Density and/or Precedence Density plots with different color modes are selected, the Color Mode field is empty (blank).
- You can select a different color mode from the dropdown menu. The selected mode is applied to all selected Density and/or Precedence Density plots.

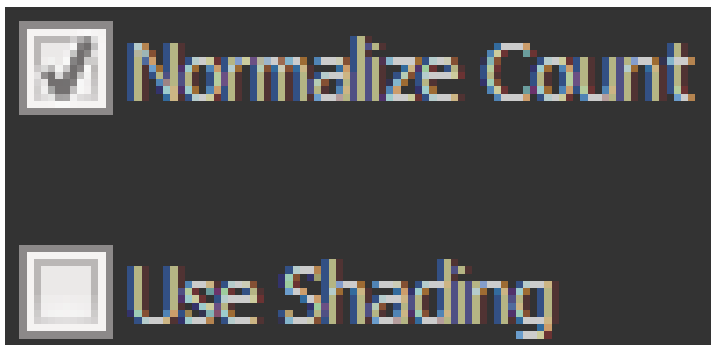
---

**Note:** The *linear mode* bins data by assigning a color index for each density pixel linearly such that each increment is determined by dividing the range ( $Z_{\max} - Z_{\min}$ ) by the number of color steps. The *log mode* bins data by assigning a color index for each density pixel logarithmically such that each increment is determined by dividing the logarithmic range ( $\log Z_{\max} - \log Z_{\min}$ ) by the number of color steps. The index is then determined as  $\text{int}((\log Z_{\text{val}} - \log Z_{\min}) / \text{increment})$ .

---

## Normalize count

**Normalize Count** checkbox controls the scaling for the selected Histogram plots.



- Normalize Count option applies only to Histogram plots and it is displayed when one or more Histogram plots are selected.
- When the checkbox is unchecked, the scale label is **Count**.
- When the checkbox is checked, the Y-axis scaling is changed to scale the plot to the Histogram peak value.  
 The scale displayed is a percentage scale from 0 to 100%, where 100% is the Histogram peak value.  
 The scale label is **Percent of Max™**.

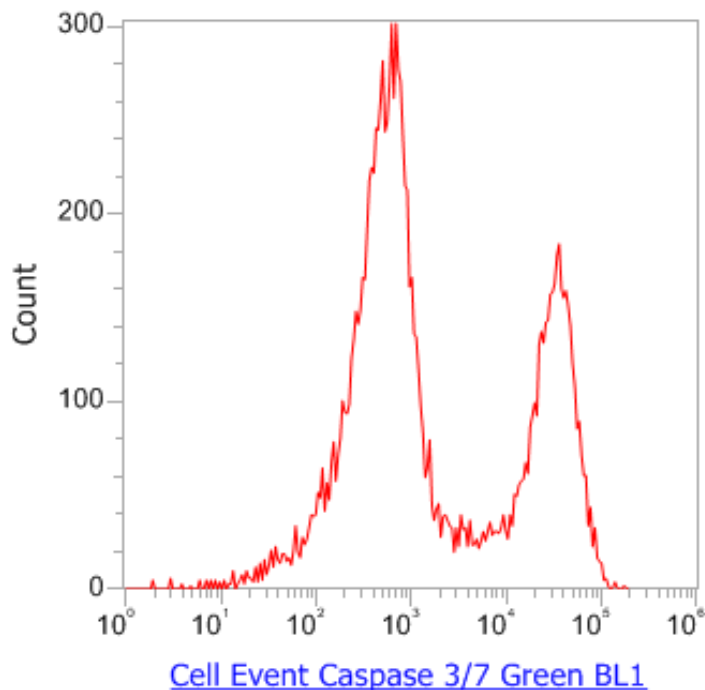


Figure 83 Normalize Count unchecked

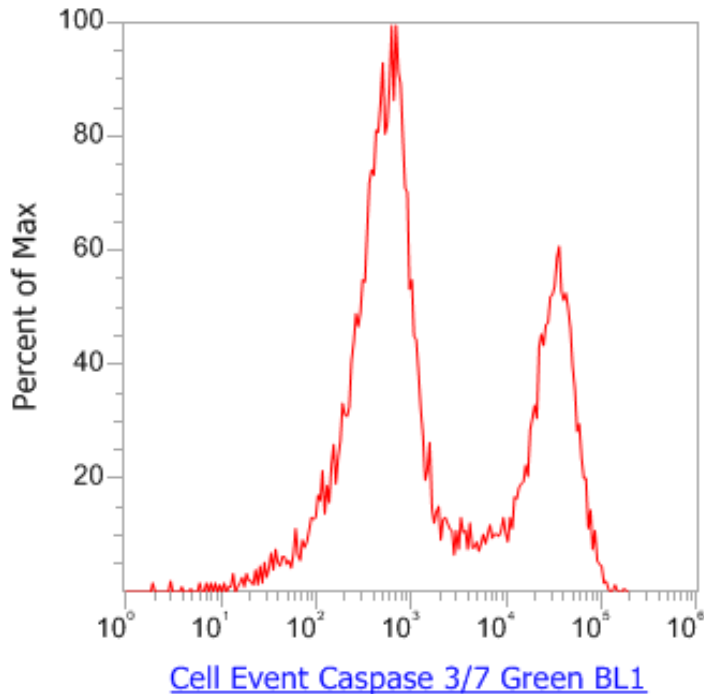
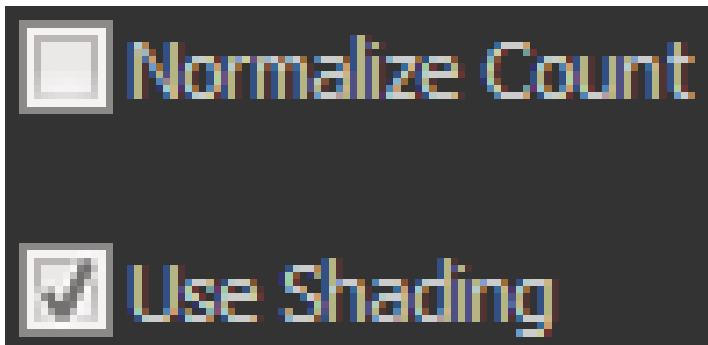


Figure 84 Normalize Count checked

### Use shading

**Use Shading** checkbox controls the shading for the selected Histogram plots.



- Use Shading option applies only to Histogram plots and it is displayed when one or more Histogram plots are selected.
- The color used for shading is the same as that used for the line color (at 64% opacity). The color and the opacity cannot be adjusted.
- When the checkbox is unchecked, shading is removed.
- When the checkbox is checked, shading is added.

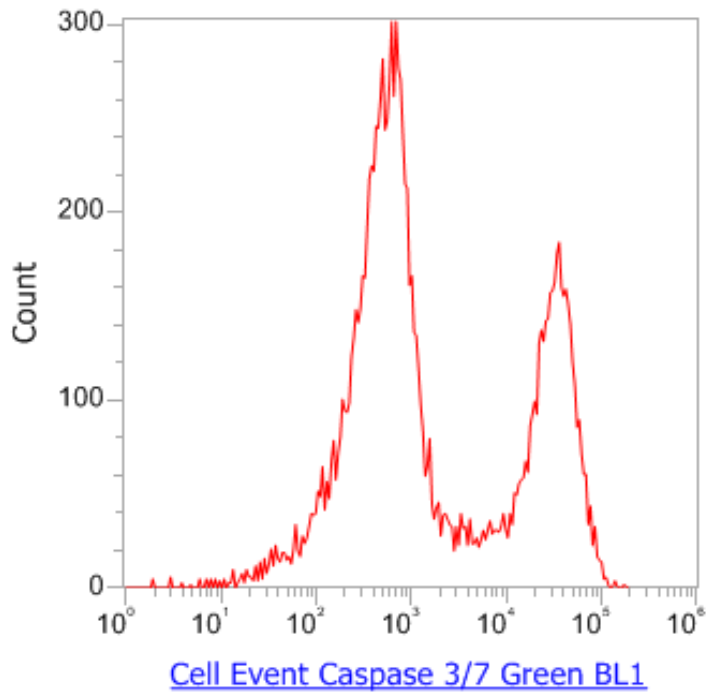


Figure 85 Use Shading unchecked

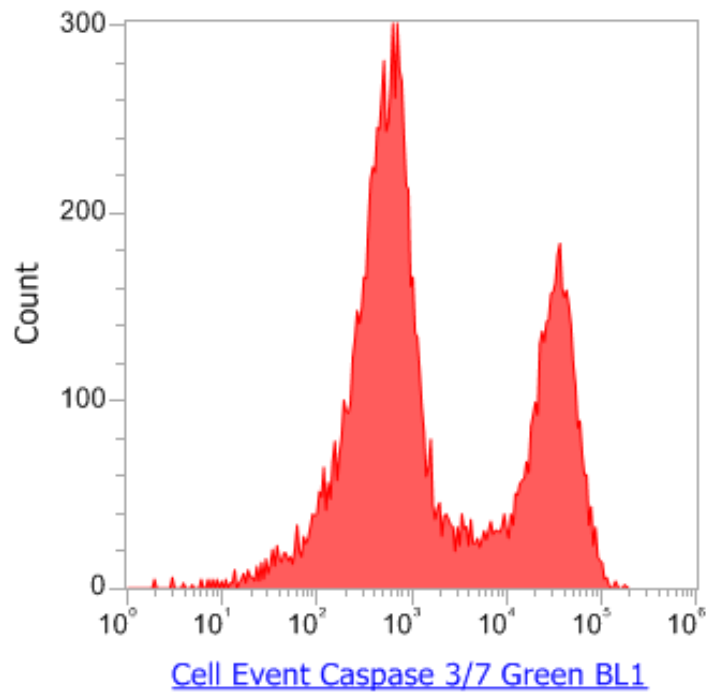


Figure 86 Use Shading checked

## Line width

**Line Width** menu controls the line thickness for the selected Histogram plots.

- Line Width option applies only to Histogram plots and it is displayed when one or more Histogram plots are selected.
- You can select or enter a number from **1 to 5**.
- If you enter an out-of-range number, the software corrects it to the nearest valid number. Non-numeric characters are not permitted.

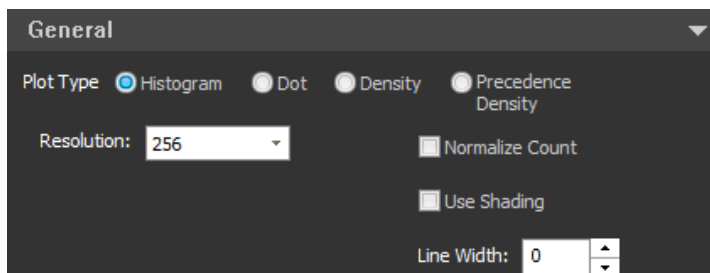
## General group – plot-specific options

This section illustrates the customization options available in the General group when single or multiple plot types are selected.

Each field in the panel displays the current settings for the selected plot. If multiple plots with differing values are selected, the value fields show the indeterminate state.

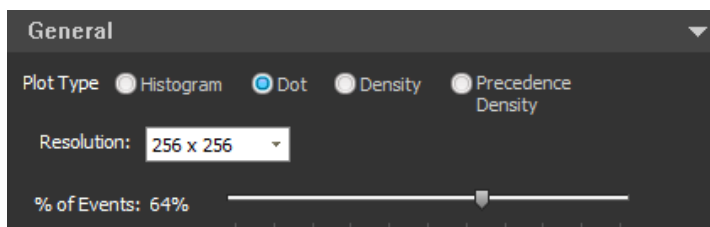
### Histogram only

- Plot Type (“Plot type radio buttons” on page 427)
- Resolution (“Resolution dropdown menu” on page 428)
- Normalize Count (“Normalize count” on page 431)
- Use Shading (“Use shading” on page 432)
- Line Width (“Line width” on page 434)



### Dot only

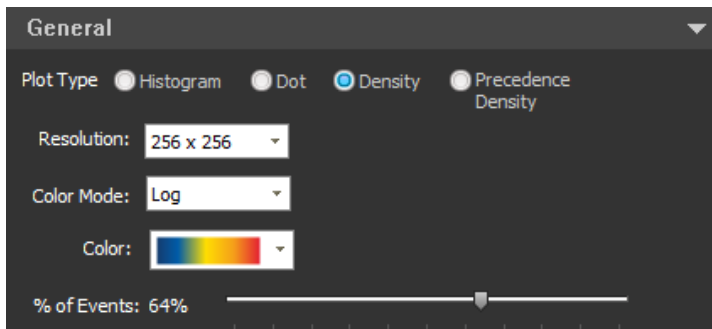
- Plot Type (“Plot type radio buttons” on page 427)
- Resolution (“Resolution dropdown menu” on page 428)
- % of Events (“% of events slider” on page 430)





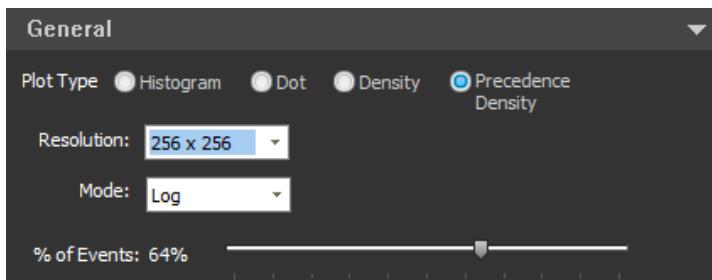
## Density only

- Plot Type (“Plot type radio buttons” on page 427)
- Resolution (“Resolution dropdown menu” on page 428)
- Mode (“Color mode dropdown menu” on page 430)
- Color (“Color dropdown menu” on page 429)
- % of Events (“% of events slider” on page 430)



## Precedence density only

- Plot Type (“Plot type radio buttons” on page 427)
- Resolution (“Resolution dropdown menu” on page 428)
- Mode (“Color mode dropdown menu” on page 430)
- % of Events (“% of events slider” on page 430)



## X-axis group

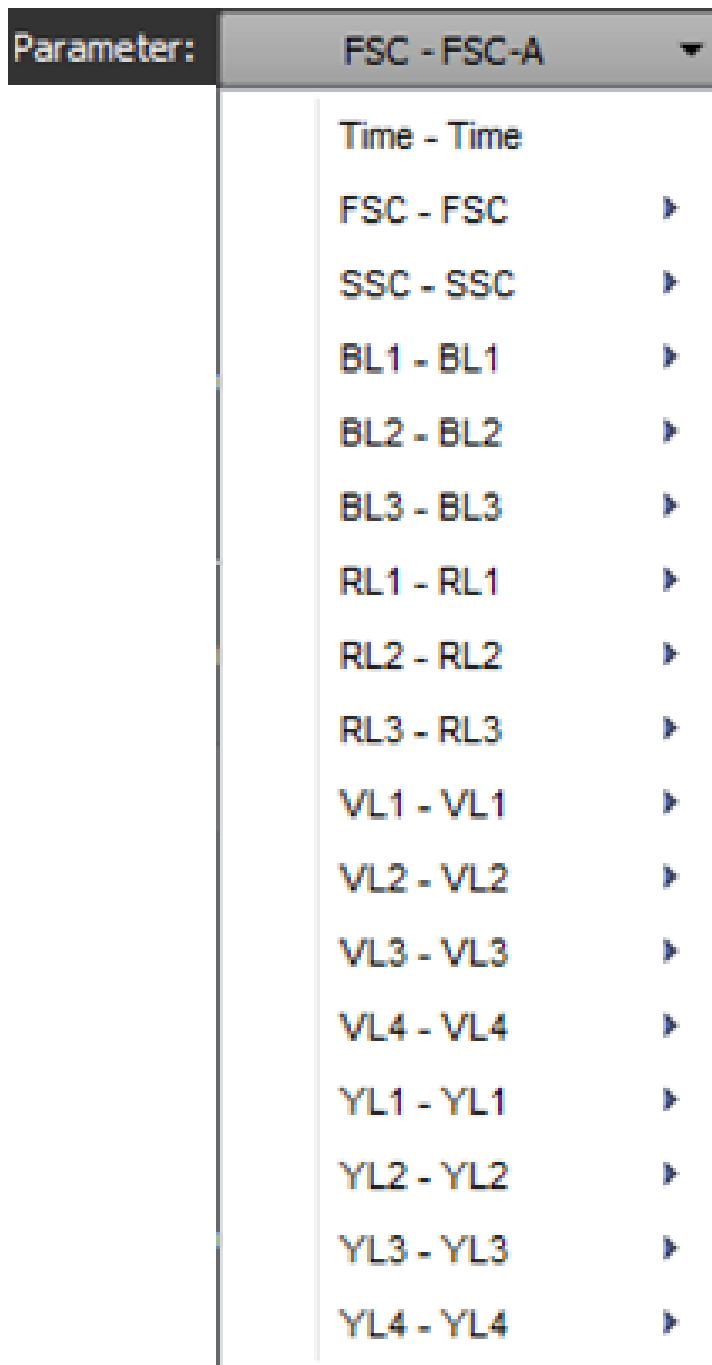
The *X-axis group* allows you to customize the X axes of selected plots.

The screenshot shows the 'X axis' customization panel. It features a title bar with 'X axis' and a dropdown arrow. Below the title bar, there are three main sections: 'Parameter', 'Scale', and 'Range'. The 'Parameter' section has a dropdown menu currently set to 'FSC-A'. The 'Scale' section has three radio button options: 'Linear' (which is selected), 'Logarithmic', and 'HyperLog™'. The 'Range' section has two radio button options: 'Automatic' and 'Manual' (which is selected). To the right of the 'Manual' option are two input fields: 'Min' with the value '1' and 'Max' with the value '1048576'.

- All customization options in the X-axis group apply to all plot types and are displayed when any plot is selected.
- The options that appear in the X-axis group vary depending on the scale selected (“Scale options” on page 439).

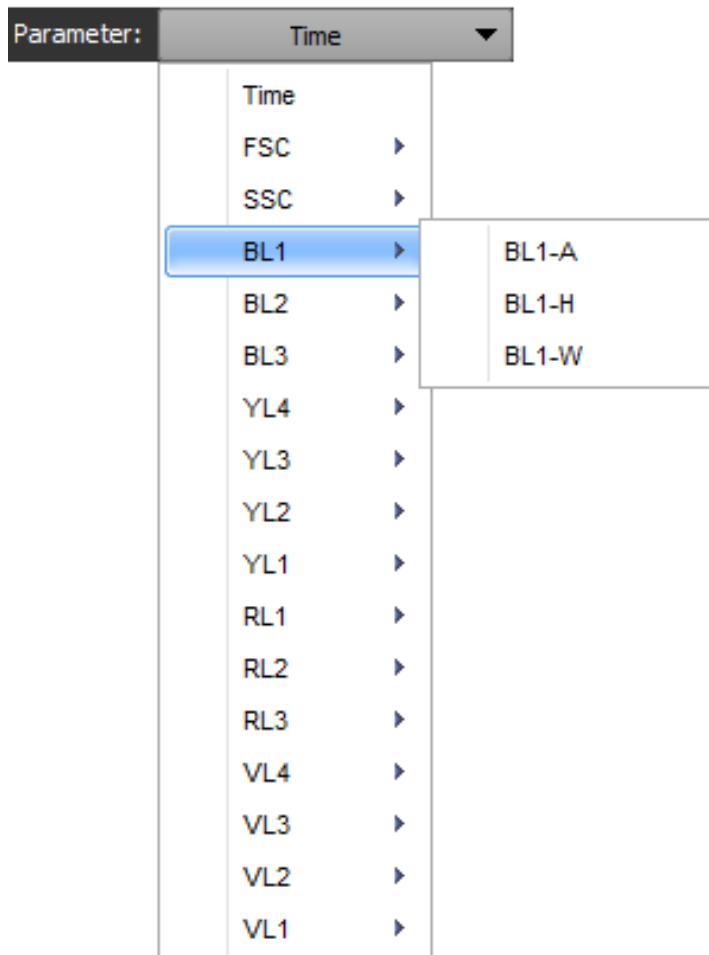
## Parameter dropdown menu

**Parameter** dropdown menu allows you to select any enabled parameter for the X-axis. All parameters in the FCS file.



- If a single plot is selected, or if multiple plots with the same X-axis parameter (including a mixture of plot types) are selected, the Parameter dropdown menu displays the parameters of the currently selected plots.

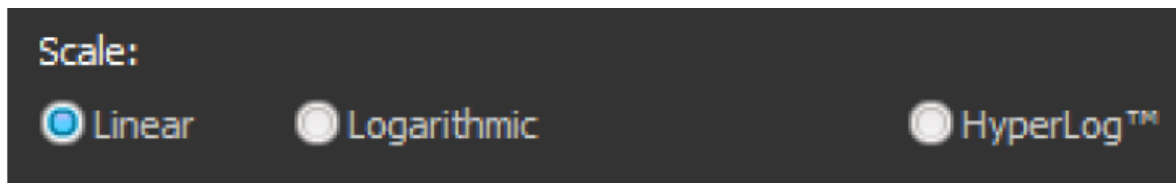
- If multiple plots of any type with different X-axis parameters are selected, the Parameter field is empty (blank).
- If the current Workspace is a Compensation Workspace, the Parameter dropdown menu is not available for dual-parameter or Histogram plots.
- To enable the Parameter dropdown menu, you can change the parameters of the gating plot within a Compensation Workspace.
- If more than 18 parameters are selected in the current Sample and if two or more measurements are enabled for the parameter, the dropdown menu includes a submenu of parameter types. Calculated parameters are always displayed in the primary menu.



- Selected plots are updated to display the selected parameter on the X-axis.

## Scale options

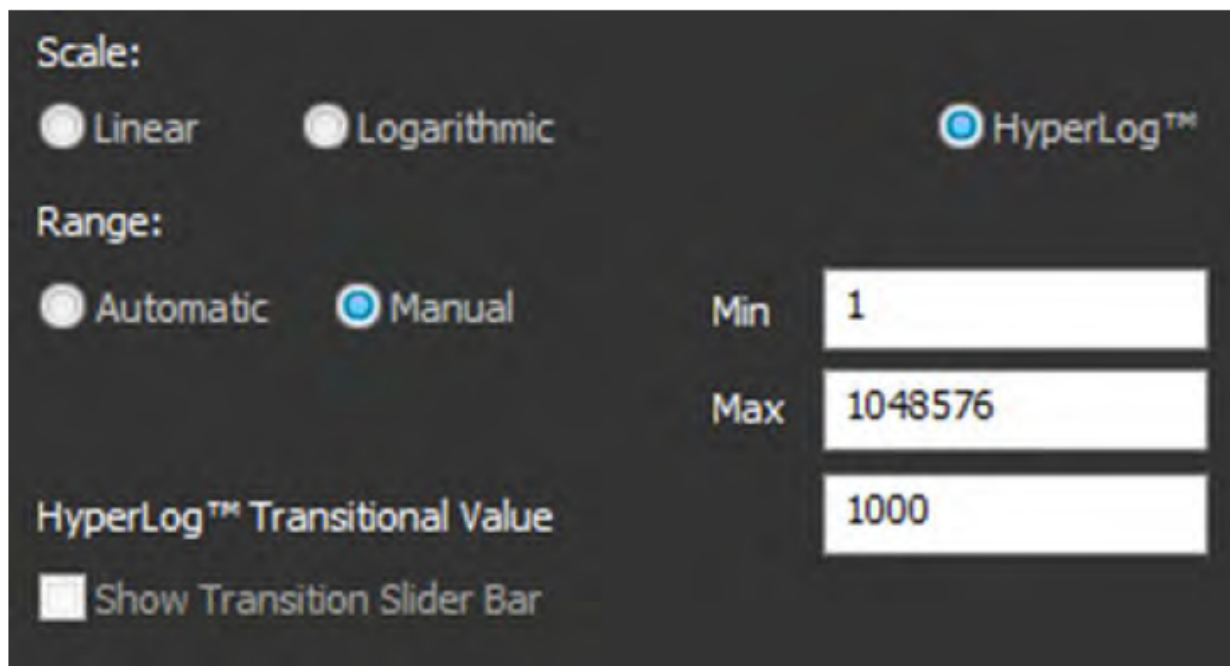
- **Scale** options allow you to select the X-axis scaling method for the selected plots: **Linear**, **Logarithmic**, and **Hyperlog™**.



**Note:** The Hyperlog™ scale uses log-linear hybrid transformations to display compensated flow cytometry data that frequently contain negative values due to compensation. Logarithmic transformations cannot properly handle negative values, and poorly display normally distributed cell types.

## Hyperlog™ scale

Hyperlog™ is an alternative bi-exponential scaling. When the **Hyperlog™** scale is selected, an additional **Transitional Value** control becomes available.



- To adjust the width of the linear area of the plot, enter a numeric value in the **Hyperlog™ Transitional Value** field. Enter a value from **0** to **1,000,000**; decimal places are not permitted.
- The Transitional Value is applied to the selected parameter.
- The Transitional Value changed on one plot changes the Transitional Value on another plot with the same parameter that is also set to Hyperlog™ scale.

## Range options

**Range** options allow you to adjust the scale range of a plot's axis.

Range:

Automatic  Manual

Min 1

Max 1048576

- **Automatic:** Sets the minimum and maximum values automatically and adjusts the upper value to the highest channel with data present.
  - For axes set to Log scaling, the minimum value is set at 1. For axes in Linear or Hyperlog™ scaling, the minimum value is set to the lowest value in the data set.
  - Data ranges are limited by the parameter range set in the FCS file (\$PnR) (see Note on page 440 below).
- **Manual:** Allows you can enter the minimum and maximum values manually. The allowable range depends on the type of scale selected.
  - **For Linear and Hyperlog™ scaling:** The default manual range is  $\pm$  the maximum range set in the FCS file (\$PnR) (see Note below).  
The allowable minimum range is  $-2^{31}$  to  $2^{31}-2$ .  
The allowable maximum range is  $-2^{31}+1$  to  $2^{31}-1$ .
  - **For Log scaling:** The default manual range is 1 to the maximum value set in the FCS file (\$PnR) (see Note below).  
The allowable minimum range is 1 to  $2^{31}-2$ .  
The allowable maximum range is 2 to  $2^{31}-1$ .
  - If **Time** is selected from the **Parameter** dropdown menu (“Parameter dropdown menu” on page 437), the Range fields are based on milliseconds.
  - If you enter an out-of-range number in the Minimum or Maximum fields, the software adjusts it to the minimum or maximum value, respectively. Non-numeric characters are not permitted.

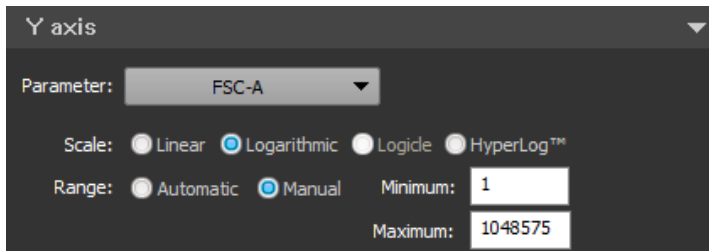
---

**Note:** \$PnR is the range for the selected parameter n (“Required keywords” on page 828). By default, this is set to  $2^{26}$  for Event count and Time parameters,  $2^{20}$  for all Height and Area measurements, and  $2^{10}$  for all Width measurements.

---

## Y-axis group

The *Y-axis group* allows you to customize the Y-axes of selected plots.



- All customization options in the Y-axis group apply to all plot types and are displayed when any plot is selected.
- The options that appear in the Y-axis group vary depending on the scale selected (see Scale options, below).

### Parameter dropdown menu

**Parameter** dropdown menu allows you to select any enabled parameter for the Y-axis. All parameters in the FCS file, including any calculated parameters, are available.

The Parameter dropdown menu for the Y-axis behaves the same as the dropdown menu for the X-axis, except when a Histogram plot is selected.

- If a Histogram plot is selected, the Y-axis parameter is fixed as **Count** and cannot be changed unless the plot is changed to a dual-parameter plot type.
- If a Histogram plot is selected and the Normalize Count checkbox is checked, the Y-axis parameter is fixed as **Percent of Max™**.
- If Histograms plots are selected along with dual-parameter plots, the Parameter dropdown menu selection has no effect on the Histogram plots.

### Scale options

**Scale** options allow you to select the Y-axis scaling method for the selected plots: **Linear**, **Logarithmic**, and **Hyperlog™**.

- **Scale** radio buttons for the Y-axis behave the same as the options for the X-axis, except when a Histogram plot is selected:
  - If a Histogram plot is selected, the Hyperlog™ scaling option is disabled.
  - If Histograms plots are selected along with two-parameter plots, all scale options are enabled; any changes to the Hyperlog™ scaling option has no effect on the Histogram plots.

### Range options

**Range** options allow you to adjust the scale range of a plot's axis. **Range** options for the Y-axis behave the same as the options for the X-axis, except when a Histogram plot is selected:

- If a Histogram plot is selected, the Y-axis is set to **Count**, the minimum value can be 0 to  $2^{31}-2$ , and the maximum value can be 1 to  $2^{31}-1$ .
- For **Percent of Max™**, the minimum value can be 1 to 99, and the maximum value can be 2 to 100.

## Text group (plots only)

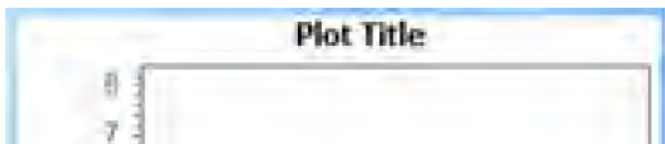
The *Text group* allows you to customize the plot text boxes. All customization options in this group apply to all plot types and are displayed when any plot is selected.

The options in the plot Text group are disabled when items other than plots are selected in the Workspace or a Compensation Workspace is displayed.

The screenshot shows the 'Text' customization panel. It includes a 'Plot Title' text field containing 'plot title', a font size dropdown set to '10', and buttons for Bold (B), Italic (I), and Underline (U). Below these are checkboxes for 'Include name in plot title' with options for Parent Gate (checked), Experiment, Group, and Sample (checked). There are also fields for 'X Axis Title' and 'Y Axis Title', another font size dropdown set to '10', and buttons for Bold, Italic, and Underline. At the bottom, there are checkboxes for 'Include name in axis title' with options for Channel Name, Target Name, and Label Name, and a 'Font' field set to 'Tahoma' with a font style dropdown set to 'AaBbC' and a font color icon.

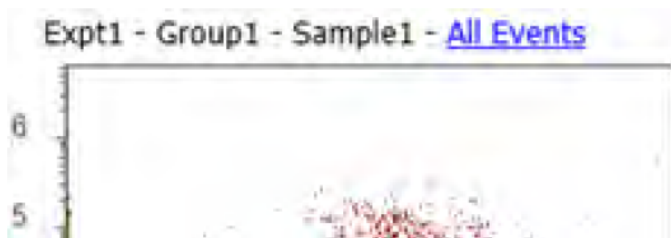
### Plot title options

**Plot title** options enable you to enter text and select font formats for the plot title. The formatted text appears at the top of all selected plots.



- You can enter up to 50 characters in the **Plot title** field; any character is permitted. If the plot title is too large to fit above the selected plots, the software truncates the title.
- In the **font size** field, you can select a font size from the dropdown or enter a number from **6** to **72**.
- You can select the **bold**, **italics**, and/or **underline** options to format the font.
- You can include the **Experiment name**, **Group name**, **Sample name**, and/or **Parent gate** by checking the appropriate **Include** checkboxes. The software appends the selections to the plot title, separated by hyphens.





## X-axis and Y-axis title options

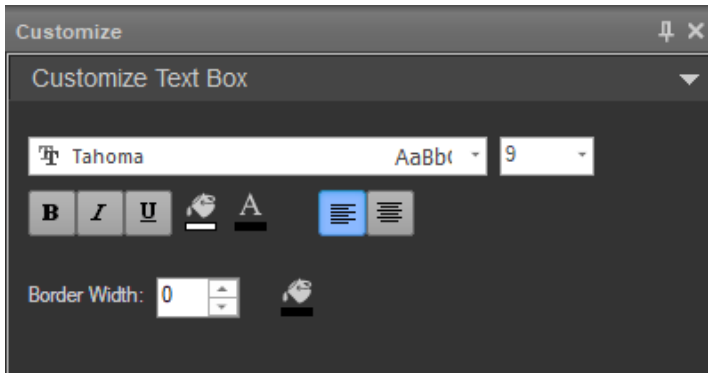
The **X-axis title** and **Y-axis title** options allow you to enter text and select font formats for the X- and Y-axis titles. The formatted text appears as the X- and Y-axis labels for all selected plots.

- You can enter up to 50 characters in the X-axis and Y-axis title fields; any character is permitted. If an axis title is too large to fit in the selected plots, the software truncates the title.
- For the X-axis:
  - The tick channel labels change from displaying the full channel number to thousands or millions, as appropriate.
  - The tick channel labels show exponential numbers if the channel number entered is >10,000.
  - The size of the labels cannot be changed.
  - Parameter names are appended with (10<sup>3</sup>) or (10<sup>6</sup>) to indicate channel scaling.
- For the Y-axis:
  - If **Histogram** plots are selected, the Y-axis title defaults to **Count** if all Histogram plots are the same and none are normalized.  
If **Histogram** plots are all set to **Normalize Count**, the title defaults to **Percent of Max**.
  - The tick channel labels show exponential numbers if the channel number entered is >10,000.
  - The size of the labels cannot be changed.
  - Parameter names are appended with (10<sup>n</sup>) to indicate channel scaling.
- In the **X-axis font size** and **Y-axis font size** fields, you can select a **font size** from the dropdown or enter a number from **6** to **72**.
- You can select the **bold**, **italics**, and/or **underline** to format the font.
- You can include the **wavelength**, **channel name**, **target name**, and/or **fluorophore name** by checking the appropriate **Include** checkboxes. The software appends the selections to the axis labels, separated by hyphens.
  - If you do not check any Include checkboxes, the axis label defaults to the parameter name specified in the **Display Parameter Name As** setting in the **General** section of the **Options dialog** (“Display Parameter Name as” on page 645).  
For example, if \$PnS (Stain Name) is selected, axis label defaults the parameter name to the stain name (for example, CD4-FITC).
  - If the target and fluorophore names are selected but unavailable (i.e., the names are from third party support/imported FCS files), the default parameter or channel name included in the imported FCS file are used.
  - Include options are not applied to the Y-axis of Histogram plots when the name is set to **Count** or **Percent of Max**, as described above.

# Customize text box options

## Text box options

*Customize Text box* group allows you to customize the font style and borders of text boxes. These options are displayed when one or more text boxes are selected in a Workspace.



## Style options

**Style** options allow you to select the font type, size, and formatting for all selected text boxes.

- You can select a **font type** from the dropdown menu. The dropdown menu includes all fonts installed on the system. The default font is Tahoma.
- You can select a **font size** from the dropdown menu or enter a number from **6** to **72**.
- You can select the **bold**, **italics**, and/or **underline** options to format the font.
- You can select to **left-align** or **center** the text.
- You can select **colors** for the text and for the Text box fill. The software uses the standard color picker dialog.

## Border options

**Border** options allow you to select the border width and color for all selected text boxes.

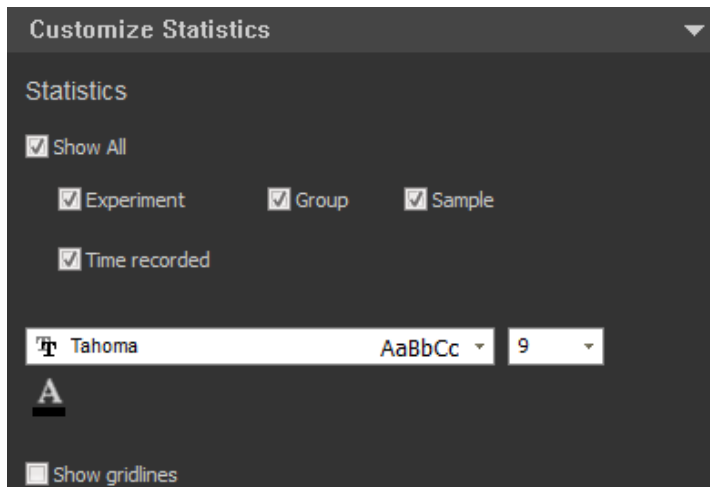
- You can select a **border width** or enter an integer from **1** to **5**.
- You can select a **color** for the border. The software uses the standard color picker dialog.

# Customize statistics box options

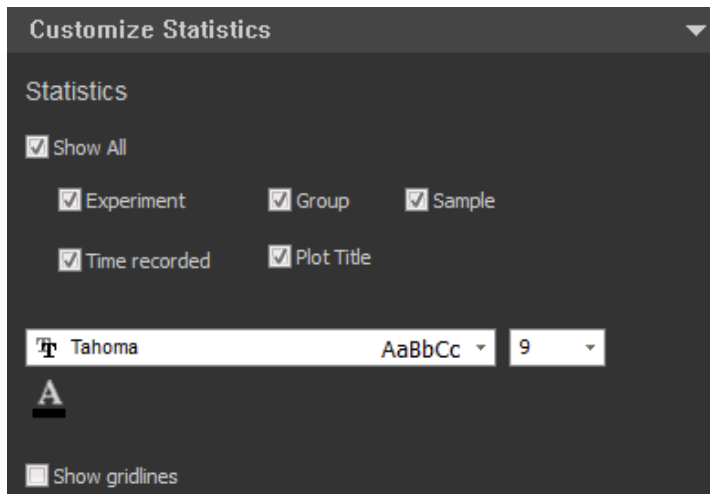
## Statistics box options

*Customize Statistics box* group allows you to customize the formatting of statistics boxes and the statistics they displayed.

- The customization options are displayed when one or more Sample-level or Plot-level statistics boxes are selected in a Workspace. The options vary slightly between sample- and Plot-level statistics boxes.
- The following option is displayed when one or more **Sample-level** statistics boxes are selected:



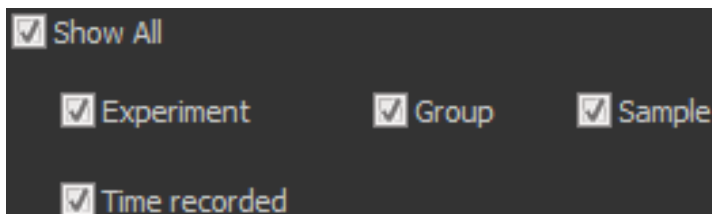
- The following option is displayed when one or more **Plot-level** statistics boxes are selected:



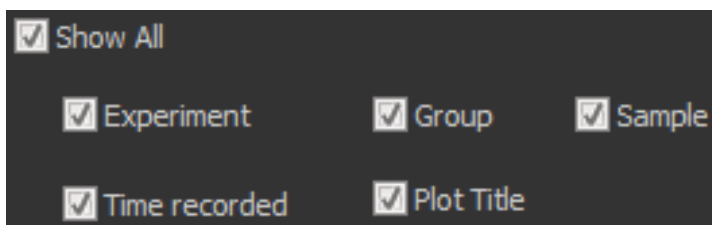
## Header options

**Header** options allow you to select which elements are automatically added to the headers of all selected statistic boxes.

- For Sample-level statistics boxes, you can include the **Experiment**, **Specimen**, and **Sample names** and the **Time Recorded** by checking the appropriate checkboxes.



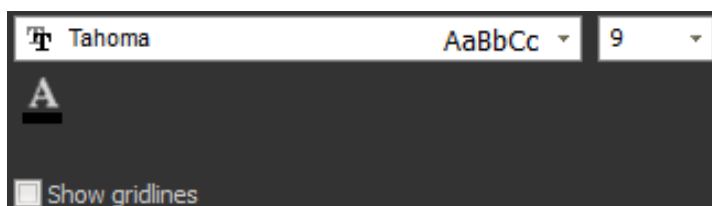
- For Plot-level statistics boxes, you can include the **Experiment**, **Group**, and **Sample** names, the **Time Recorded**, and the **Plot title** by checking the appropriate checkboxes.



- Checking the **Show All** checkbox selects all available elements.
- The software includes your selections in the statistics box header.

## Formatting options

**Formatting** options allow you to select the font type, size, and formatting for all selected statistics boxes.



- You can select a **font type** from the dropdown menu. The dropdown menu includes all fonts installed on the system. The default font is Tahoma.
- You can select a **font size** from the dropdown menu or enter a number from **6** to **72**.
- You can select the **bold**, **italics**, and/or **underline** options to format the font.
- You can select a **color** for the text. The software uses the standard color picker dialog.
- You can display a border around the box and gridlines inside the box by checking the **Show Gridlines** checkbox. The border and gridline colors and widths cannot be adjusted.

# Customize gate options

## Overview

*Customize Gate options* allow you to customize the properties of the following gate types:

- Rectangle, Oval, and Polygon
- Contour Autogate and Ellipse Autogate
- Histogram
- Quadrant

---

**Note:** To set default options for gates, use the *Gate Options dialog* (“Gate Options” on page 656).

---

## Customize panel behavior for gates

- The Customize Gate options are displayed only when a gate is selected.
- The Customize Gate options are arranged into the following groups:
  - *Gate Type* (“Gate type group” on page 448)
  - *Name options* (“Name options” on page 459)
  - *Coordinates* (“Coordinates group” on page 461)
  - *Autogating* (available only for Contour and Ellipse Autogates) (“Autogating options” on page 464)
- Each group is contained in a group box that is collapsible.
- The options displayed in each group vary, depending on the gate types and options selected.
- The values displayed in any field are the current settings for the selected gate.
- When a Compensation sample is active, all gate customization options are disabled, except for the options specified in Chapter 21, “Compensation”.

## Gate type group

*Gate Type group* controls the gate type and general gate characteristics (such as opacity and color). The customization options that appear in the Gate Type group vary, depending on the gate types selected.

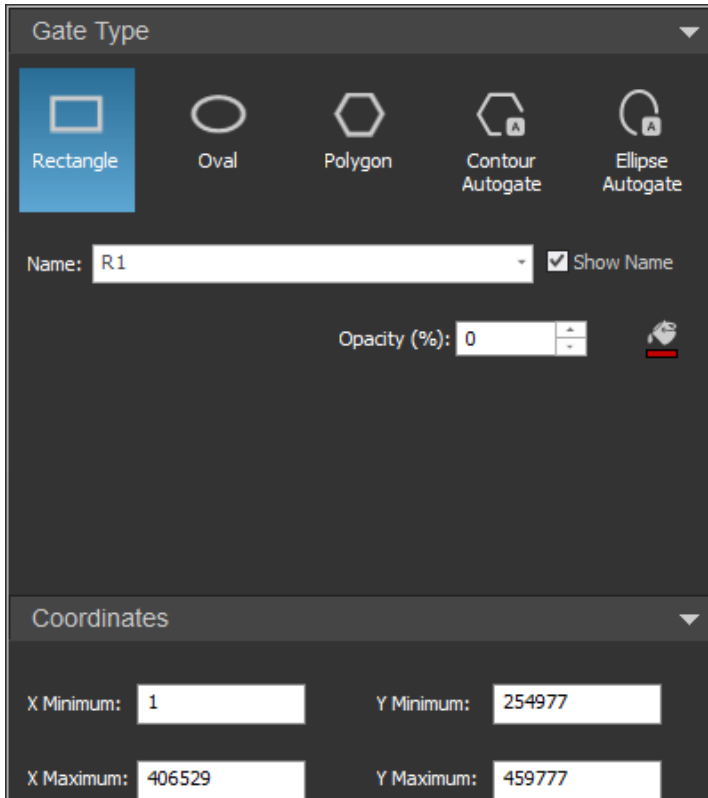


Figure 87 Rectangle gates

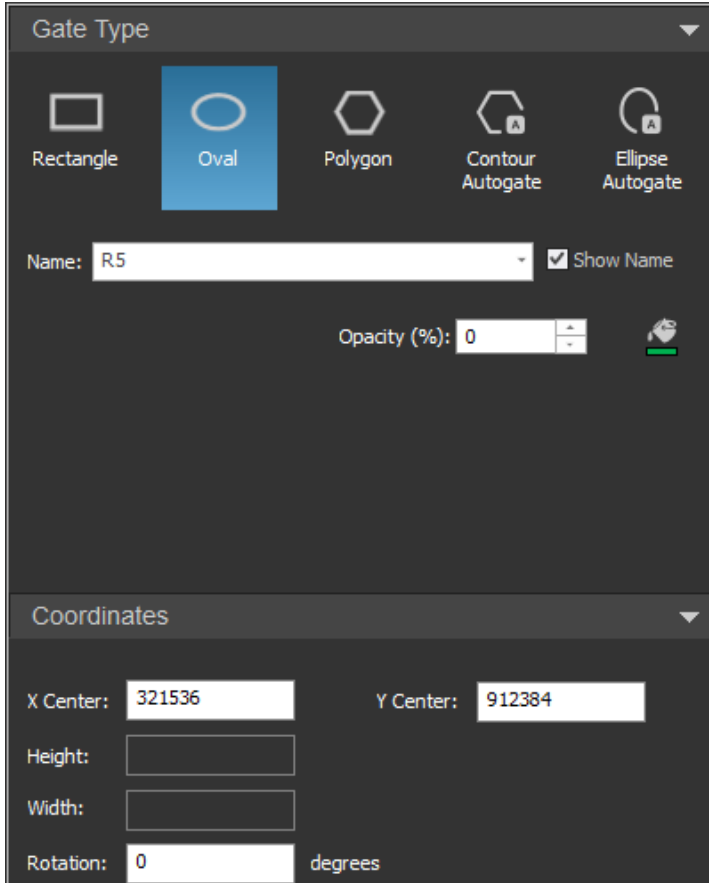


Figure 88 Oval gates

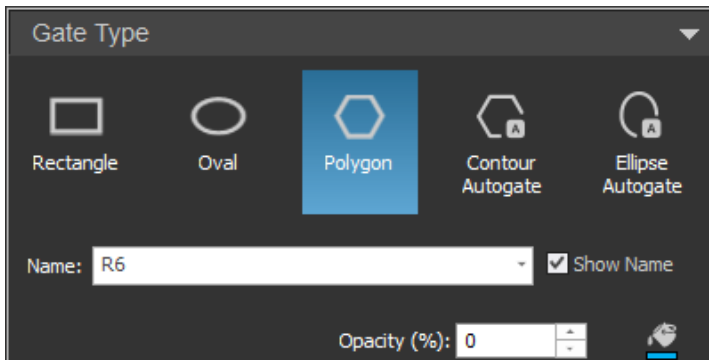


Figure 89 Polygon gates

The image shows a software interface for configuring a 'Contour Autogate'. At the top, there is a 'Gate Type' dropdown menu with a blue arrow pointing down. Below it are five icons representing different gate shapes: Rectangle, Oval, Polygon, Contour Autogate (which is highlighted with a blue background), and Ellipse Autogate. Below the icons is a 'Name' field containing 'R.2' and a 'Show Name' checkbox which is checked. There are four main configuration sections: 1) 'Autogating on' checkbox (checked) and 'Opacity (%)' field (0); 2) 'Preset' dropdown (3) and 'Min events' field (200); 3) 'Density level' field (6) and 'Resolution' field (128); 4) A 'Reset' button. At the bottom, there is a 'Coordinates' dropdown menu with a blue arrow pointing down. Below it are four input fields: 'X Minimum' (422913), 'Y Minimum' (403457), 'X Maximum' (832512), and 'Y Maximum' (608256).

Figure 90 Contour Autogate



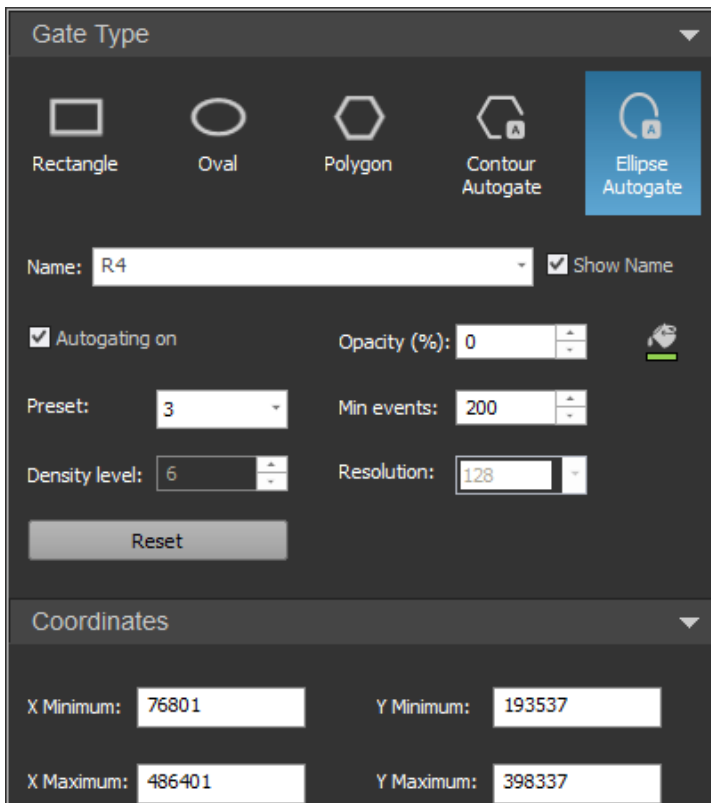


Figure 91 Ellipse Autogate

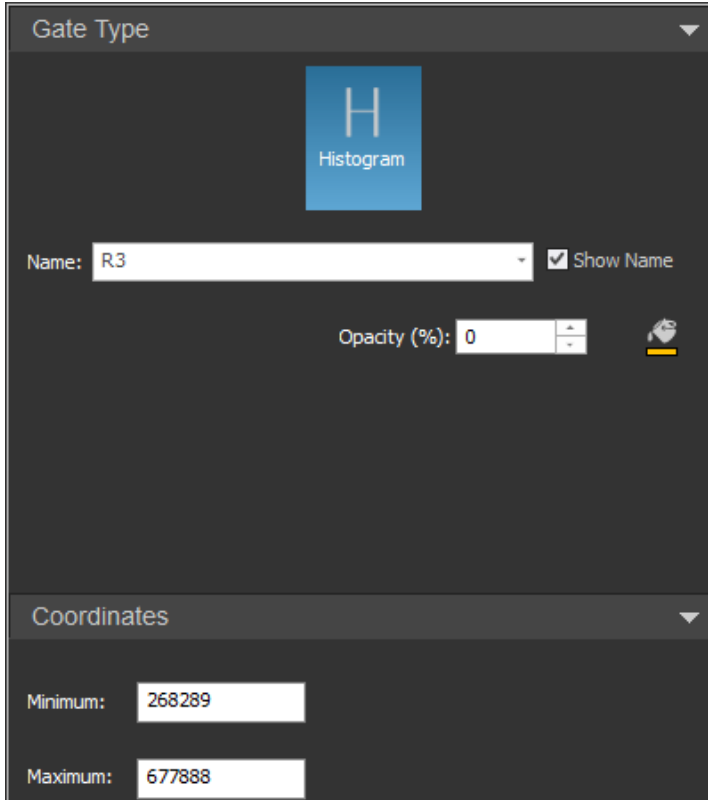


Figure 92 Histogram gates

Gate Type

Quad Bent Quad

Use smart gate naming

UL Name   Show name

Opacity %

UR Name   Show name

Opacity %

LL Name   Show name

Opacity %

LR Name   Show name

Opacity %

Coordinates

X Center  Y Center

Link Gates

Gate

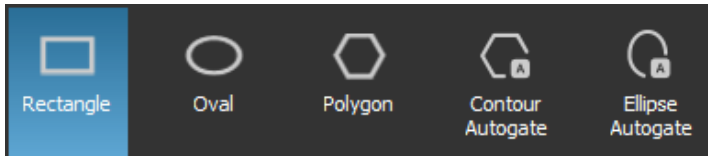
Figure 93 Quadrant gates

## Gate type buttons

The Gate Type buttons allow you to change the gate type of the selected gate on a plot. The button for the currently selected gate type is highlighted.

### Rectangle, oval, and polygon gates

When a **Rectangle**, **Oval**, **Polygon gate**, **Contour Autogate**, or **Ellipse Autogate** is selected, all gate type buttons are displayed together. The button for the selected gate type is highlighted.



- To change the gate type of the selected gate, click the desired **Gate Type** button.
- The following figures show the same area of the plot with different gate types.

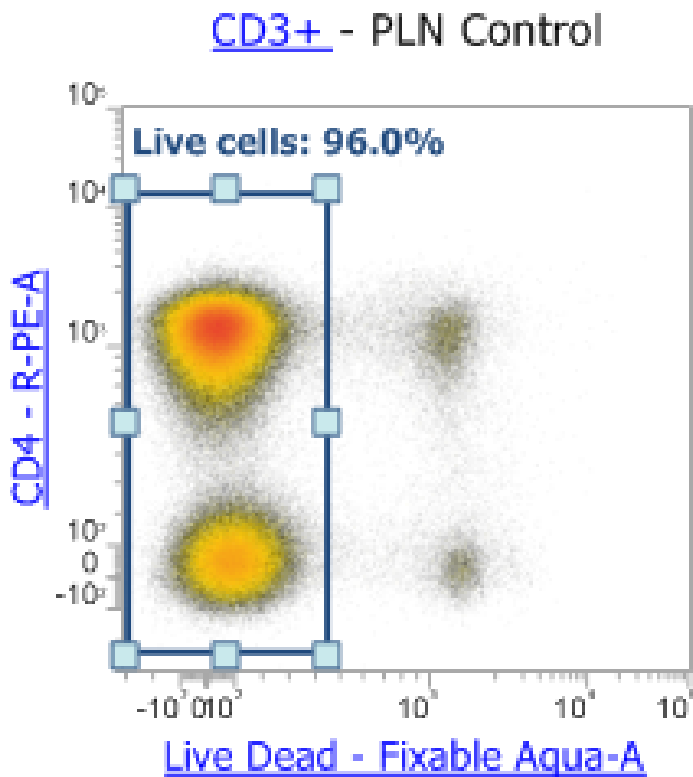


Figure 94 Rectangle gate

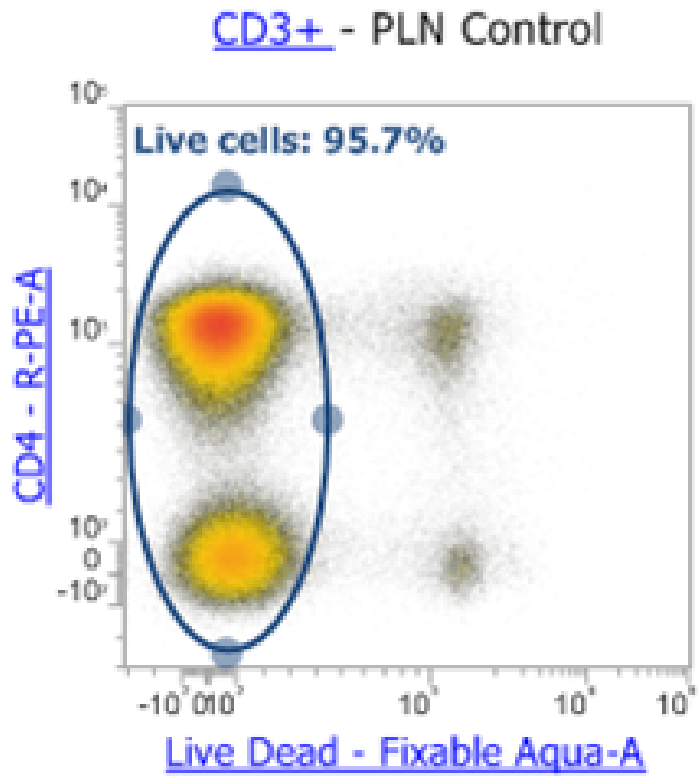


Figure 95 Oval gate

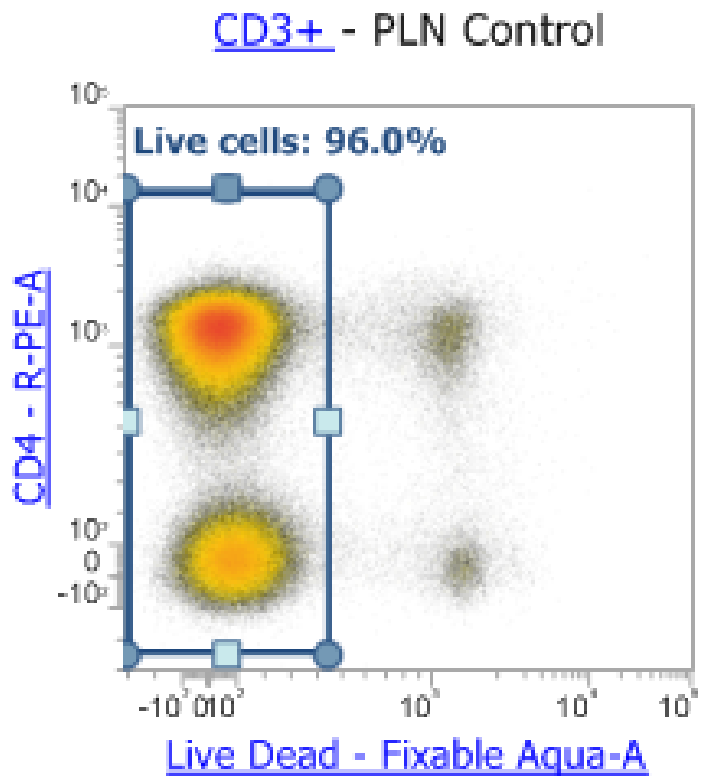
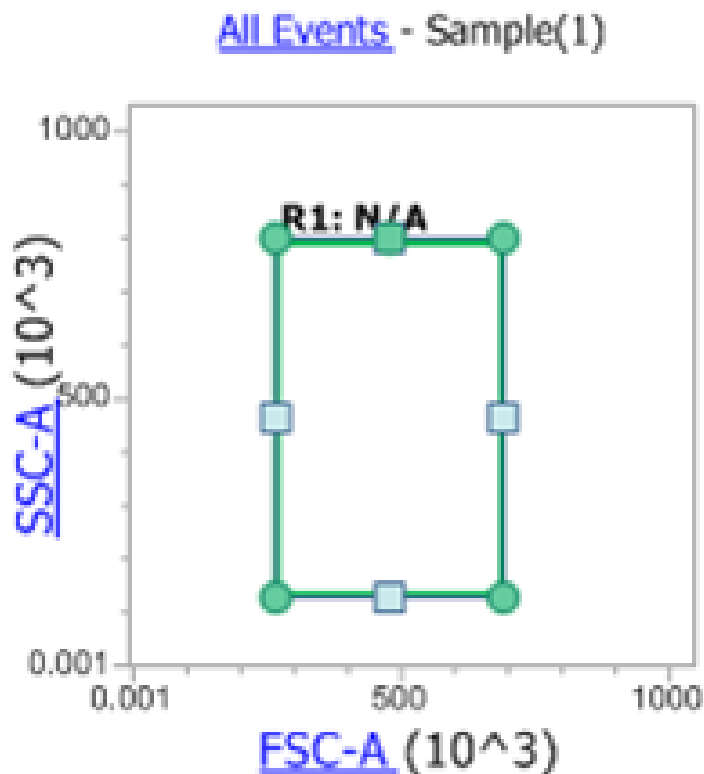


Figure 96 Polygon gate

- When you switch to a different gate type, the software:
  - Creates a new gate with the same boundaries as the original gate (X minimum, X maximum, Y minimum, and Y maximum)
  - Sizes the new gate according to the original gate's size and position.
- When you switch to a Polygon gate, the software creates a 5-point polygon in the shape of the bounding rectangle as shown in the following example.



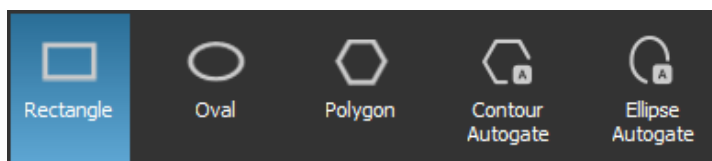

---

**Note:** For a description of the behaviors when you switch to and from a Contour Autogate or Ellipse Autogate to other gate types, see “Contour autogate and ellipse autogate” on page 456.

---

### Contour autogate and ellipse autogate

When a **Rectangle**, **Oval**, **Polygon gate**, **Contour Autogate**, or **Ellipse Autogate** is selected, all gate type buttons are displayed together. The button for the selected gate type is highlighted.



To change the gate type of the selected gate, click the desired **Gate Type** button.

The following tables describe the behavior when a Contour or Ellipse Autogate is switched to another gate type and vice versa.

**Table 1 Conversion of Contour Autogate to another gate**

New gate	Behavior
Rectangular gate	Bounding box of the autogated ellipse generates a new gate.
Oval gate	
Polygon gate	Polygon vertices persist.
Ellipse autogate	New gate is generated based on polygon vertices. Smoothing vertices and other smoothing behaviors are applied.

**Table 2 Conversion of other gates to Contour Autogate**

Old gate	Behavior
Rectangular gate	Bounding box of the Rectangle gate generates a Contour autogate bounding box.
Oval gate	Bounding box of the Oval gate generates a Contour autogate bounding box.
Polygon gate	Bounding box of the Polygon gate generates a Contour autogate bounding box.
Ellipse autogate	A Contour autogate is generated based on the Ellipse autogate bounding box and polygon vertices with same scanning region. Smoothing behavior is <b>not</b> applied.

**Table 3 Conversion of Ellipse Autogate to another gate**

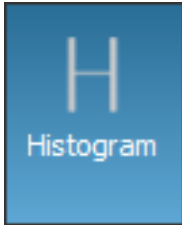
New gate	Behavior
Rectangular gate	Bounding box of the autogated polygon generates a new gate.
Oval gate	
Polygon gate	New gate is generated based on polygon vertices. Smoothing vertices and other smoothing behaviors are <b>not</b> applied.
Contour autogate	

**Table 4 Conversion of other gates to Ellipse Autogate**

Old gate	Behavior
Rectangular gate	Bounding box of the Rectangle gate generates an Ellipse autogate bounding box.
Oval gate	Bounding box of the Oval gate generates an Ellipse autogate bounding box.
Polygon gate	Bounding box of the Polygon gate generates an Ellipse autogate bounding box.
Contour autogate	An Ellipse autogate is generated based on the Contour autogate bounding box and polygon vertices with same scanning region.

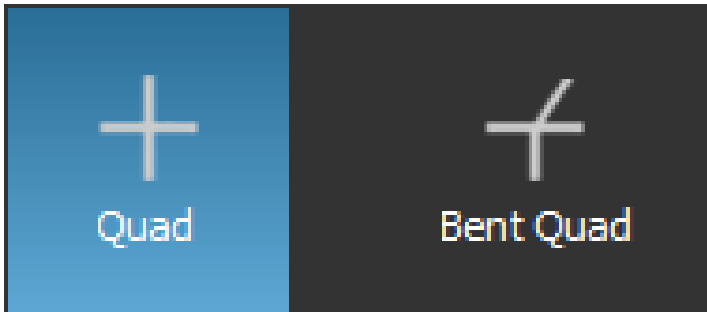
## Histogram gates

When a **Histogram gate** is selected, only the Histogram button is displayed. You cannot change a Histogram gate to a different gate type.



## Quadrant gates

When a **Quadrant gate** is selected, you can choose between the **Orthogonal** and **Bent Quadrant** gates.



- Orthogonal mode maintains the quadrant orthogonally as the center point is moved. In the Bent mode, all vertices move independently of each other.
- The Quadrant mode selection is mutually exclusive. Selecting the Orthogonal Quadrant option deselects the Bent Quadrant option and vice versa.
- If the end points of a quadrant gate are moved on the Workspace or Filmstrip view, the quadrant gate is automatically set as a Bent Quadrant gate.



## Name options

### Name field

**Name field** displays the name of the currently selected gate. The Name options allow you to assign a gate name and show or hide the gate name on the plot.



The screenshot shows a user interface for naming a gate. It consists of a text input field labeled 'Name:' containing the text 'R5'. To the right of the input field is a dropdown menu with a downward-pointing arrow, which is currently open, showing three options: 'Lymphs', 'Grans', and 'Monos'. To the right of the dropdown menu is a checkbox labeled 'Show Name', which is checked.

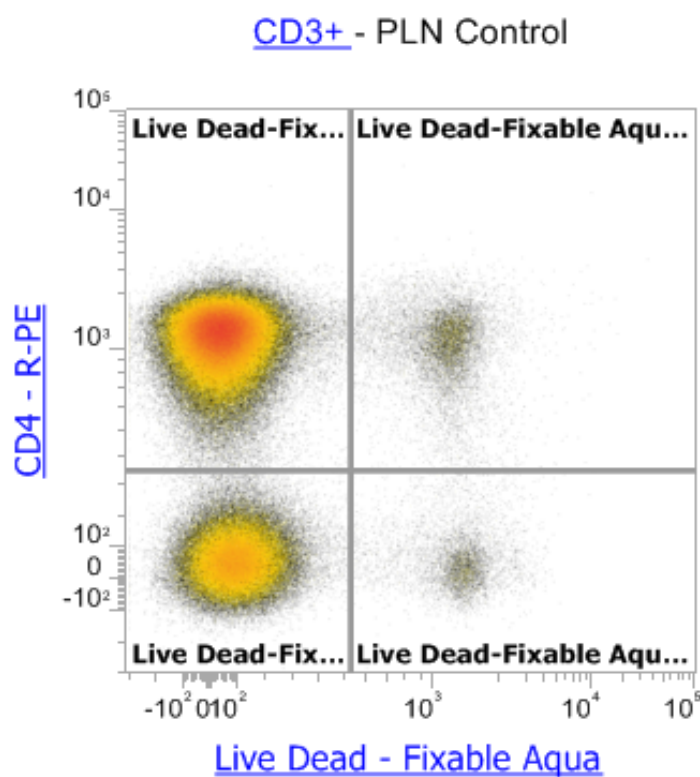
- You can enter up to 50 characters in the Name field; any character is permitted. Alternatively, you can select a gate name from the **Quick Select** dropdown menu. The dropdown menu options are set in *Gate Options* in the *Options dialog* (“Display Parameter Name as” on page 645).
- The gate name must be unique. If you enter a duplicate name, the software appends a numerical suffix in parentheses; for example, **GateName (2)**.
- You can show or hide the gate name on the plot by checking or unchecking the **Show Name** checkbox.
- For Quadrant gates, four Name fields are displayed (one for each quadrant).
- For Quadrant gates, you can select the **Use smart gate naming** option (see below) instead of entering the gate names.

## Smart gate naming

**Smart Gate Naming** option allows you to select and deselect automatic gate naming. This option is available only for Quadrant gates.



- If you check the **Use smart gate naming** check box, the software automatically names the Quadrant gates, using the options specified in *Gate Options* in the *Options dialog* (“Display Parameter Name as” on page 645).



- For Quadrant gates, the four gate names are generated (one for each quadrant).
- Smart names can be up to 50 characters; any character is permitted.
- Smart names must be unique. If a duplicate name is generated, the software appends a numerical suffix in parentheses; for example, **GateName (2)**.
- When you edit a smart name, **Use smart gate naming** is automatically unchecked. You must enter the gate names manually, as described in **Name field** above.
- When you uncheck **Use smart gate naming**, you must enter the gate names manually, as described in **Name field** above.

---

**Note:** Smart Gate Naming option is available only for Quadrant gates.

---

## Coordinates group

The customization options that appear in the Coordinates group vary depending on the gate type selected. This section describes all possible options available in the Coordinates group.

---

**Note:** Coordinate options are not available for Polygon gates.

---

### Coordinates group for rectangular gates, contour autogates, and ellipse autogates

For a Rectangular gate or the bounding box of Contour or Ellipse autogates, you can enter **X** and **Y** **minimum** and **maximum coordinates**.

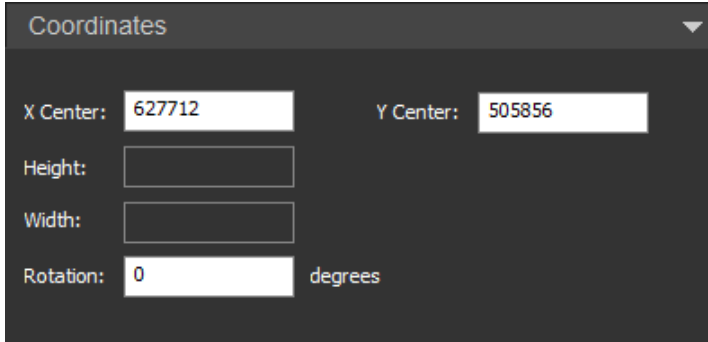


The screenshot shows a dark-themed panel titled "Coordinates" with a dropdown arrow on the right. It contains four input fields arranged in a 2x2 grid. The top row has "X Minimum:" followed by a text box containing "422913" and "Y Minimum:" followed by a text box containing "403457". The bottom row has "X Maximum:" followed by a text box containing "832512" and "Y Maximum:" followed by a text box containing "608256".

- Non-numeric characters are not permitted.
- The allowable X and Y coordinates depend on the type of scale selected.
  - **Linear and Hyperlog™ scaling:**  
The allowable minimum range is  $-2^{31}$  to  $2^{31}-2$ .  
The allowable maximum range is  $-2^{31}+1$  to  $2^{31}-1$ .
  - **Log scaling:**  
The allowable minimum range is 1 to  $2^{31}-2$ .  
The allowable maximum range is 2 to  $2^{31}-1$ .
- If you enter an out-of-range number, the software defaults to the minimum or maximum value in the allowable range.

## Size and coordinates group for oval gates

For an Oval gate, you can enter **center point coordinates**, and the **height**, **width**, and **rotation** values.



Coordinates

X Center:  Y Center:

Height:

Width:

Rotation:  degrees

- Non-numeric characters are not permitted.
- The allowable X and Y coordinates depend on the type of scale selected:
  - **Linear and Hyperlog™ scaling:** The allowable range is  $-2^{31}$  to  $2^{31}-1$ .
  - **Log scaling:** The allowable range is 1 to  $2^{31}-1$ .
- The allowable height range is  $-2^{31}$  to  $2^{31}-1$ .
- The allowable width range is  $-2^{31}$  to  $2^{31}-1$ .
- The allowable rotation range is  $-360$  to  $360$ .
- If you enter an out-of-range number, the software defaults to the minimum or maximum value in the allowable range.

## Coordinates group for histogram gates

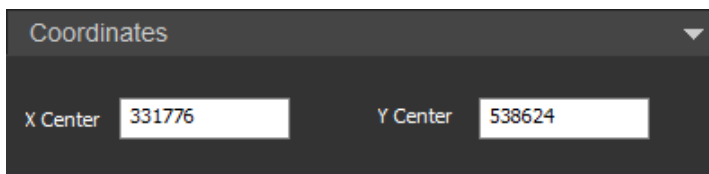
For a Histogram gate, you can enter **left** and **right boundary coordinates**.



- Non-numeric characters are not permitted.
- The allowable coordinates depend on the type of scale selected:
  - **Linear and Hyperlog™ scaling:**  
The allowable minimum range is  $-2^{31}$  to  $2^{31}-2$ .  
The allowable maximum range is  $-2^{31}+1$  to  $2^{31}-1$ .
  - **Log scaling:**  
The allowable minimum range is 1 to  $2^{31}-2$ .  
The allowable maximum range is 2 to  $2^{31}-1$ .
- If you enter an out-of-range number, the software defaults to the minimum or maximum value in the allowable range.

## Coordinates group for quadrant gates

For a Quadrant gate, you can enter the **center point coordinates**.



- Non-numeric characters are not permitted.
- The allowable X and Y coordinates depend on the type of scale selected:
  - **Linear and Hyperlog™ scaling:** The allowable range is  $-2^{31}$  to  $2^{31}-1$ .
  - **Log scaling:** The allowable range is 1 to  $2^{31}-1$ .
  - If you enter an out-of-range number, the software defaults to the minimum or maximum value in the allowable range.

## Autogating options

Autogating options in the Customize Gate Options panel allow you to customize settings for the selected Contour or Ellipse Autogate.

Customize Autogating options are available only when a Contour Autogate or an Ellipse Autogate is selected from a plot in the Workspace.

**Note:** To set default Autogating options, use the Options dialog ► Gate Options tab (“Gate Options” on page 656).

### Autogating on

**Autogating on** checkbox allows you to turn the autogating function for the selected auto region on or off. This control is enabled when an auto region is selected from a plot in the Workspace.

- **Autogating on** control sets the selected auto region to the following state, depending on whether it is checked or unchecked:

Previous state	New state when checked	New state when unchecked
Autogate ON	Autogate ON	Autogate OFF
Autogate ON – [NO FIT]	Autogate ON – [NO FIT] <sup>[1]</sup>	Autogate OFF
Autogate OFF	Autogate ON – [NO FIT]	Autogate OFF

<sup>[1]</sup> An autogate that is ON but with NO FIT becomes an ON autogate at this point, if the criteria for a successful autogate are met.

### Reset

**Reset** button is enabled when an auto region is selected. When clicked, autogates will revert to the global default settings set in the Gate Options menu.

## Preset

**Preset** dropdown allows you to set the **Preset** for individual autogates selected from a plot in the Workspace. A Preset consists of Density Level and Resolution settings.

- The Preset dropdown has the following options:
  - **1** (Density level: 30, Resolution: 128)
  - **2** (Density level: 10, Resolution: 128)
  - **3** (Density level: 6, Resolution: 128)
  - **4** (Density level: 1, Resolution: 128)
  - **Custom** (By default, Density level: 5 and Resolution: 64, but you can change the custom settings using the **Density level** and **Resolution** controls)
- When a Preset is selected for an autogate, the Density level and Resolution controls display the values that correspond to the Preset choice for that autogate.
- By default, Preset is set at 3. However, you can change the default Preset in the Options dialog ► Gate Options tab (“Gate Options” on page 656).

## Min events

Sets the **Min events** value for the selected autogate, which is the number of events that must be contained within the auto region for the population to be autogated.

- By default, Min events is set to 200.
- The range of values for Min events is 1 to 9,999,999.
- If you enter a Min events value that is outside the numeric limits, the value defaults to the maximum allowed number, if the value exceeds the upper limit, or to the minimum allowed value, if the entered value is less than the lower limit.

## Density level

Sets the default **Density level** for the selected autogate.

- The Density level field is populated the value that corresponds to the selected Preset option (see Preset, above).
- The range for Density level is 1 to 100.
- The Density level control is disabled unless the **Custom** option is selected in the Preset dropdown.

## Resolution

Sets the default **Resolution** for the selected autogate.

- The Resolution field is populated with the value that corresponds to the selected Preset option (see Preset, above).
- The Resolution dropdown has the following options: **64, 128, 256, 512, 1024**.
- The Resolution control is disabled unless the **Custom** option is selected in the Preset dropdown.

## Customize experiment (Plate or Tube) options

### Overview

*Customize Plate Experiment* and *Customize Tube Experiment* options allow you to customize the following:

- Plate information (available only for Plate Experiments)
- Experiment information
- Group information
- Sample information



## Customize panel for plate experiments

The screenshot displays a dark-themed interface with four vertically stacked, expandable sections, each with a downward-pointing arrow on the right side of its header:

- Plate Information:** Contains a 'Name' field with the value 'Plate(2)', a 'Type' dropdown menu set to '96 Well Round/U Bottom', two checkboxes for 'Use Foil Cover' and 'Use Cooling Block' (both are unchecked), an 'ID' field, and a 'Notes' field.
- Experiment Information:** Contains a 'Name' field with the value 'Plate' and a small icon to its right, and a 'Notes' field.
- Group Information:** Contains a 'Name' field and a 'Notes' field.
- Sample Information:** Contains a 'Name' field and a 'Notes' field.

- Customize Plate options are available only when a Plate Experiment is displayed and the *Heat Map view* (Chapter 6, “Heat Map View”) is in focus.
- The Customize Plate options are arranged into the following groups:
  - *Plate information* (“Plate information” on page 468)
  - *Experiment information* (“Experiment information” on page 470)
  - *Group information* (“Group information” on page 471)
  - *Sample information* (“Sample information” on page 473)
- Each group is contained in a group box that is collapsible.
- The values displayed in any field are the current settings.
- Plate information is automatically saved and applied to all Samples in the Plate, including any Tubes added to the Plate Experiment, after the focus is lost from the selected field.
- Experiment information is applied to the selected Experiment after the focus is lost from the selected field.

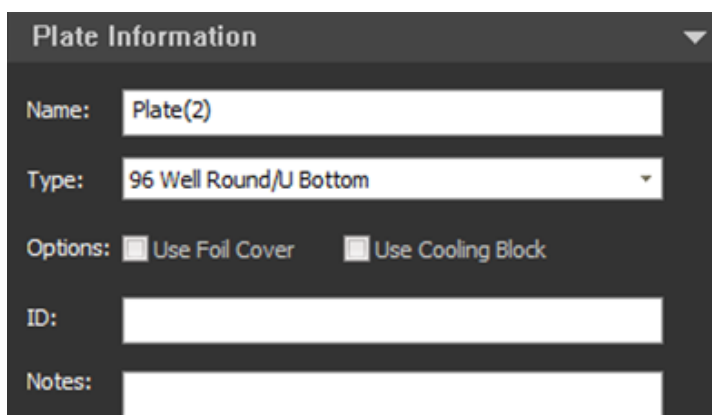
- Group information is automatically saved in the selected Group after the focus is lost from the selected field.
- Sample information is automatically saved and applied to all selected Samples after the focus is lost from the selected field.

## Customize panel behavior for tubes

When a Tube Experiment is loaded, the customizations options are the same as the options available for Plate Experiments, but the Plate information options (“Plate information” on page 468) are not displayed.

## Plate information

*Plate Information* options are available only when a Plate Experiment is displayed in the application and the Heat Map view (Chapter 6, “Heat Map View”) is in focus.



The screenshot shows a dark-themed panel titled "Plate Information" with a dropdown arrow on the right. It contains the following fields and options:

- Name:** A text input field containing "Plate(2)".
- Type:** A dropdown menu showing "96 Well Round/U Bottom".
- Options:** Two checkboxes, both of which are unchecked: "Use Foil Cover" and "Use Cooling Block".
- ID:** An empty text input field.
- Notes:** An empty text input field.

---

**Note:** Plate information options are not available when a Tube Experiment is loaded.

---

## Name

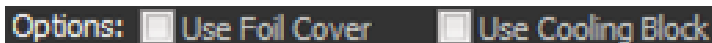
- The **Name** field displays the current name of the Plate.
- To modify the name of the selected Plate, enter the new name in the **Name** field. The entered name is automatically saved and applied to the Plate once the focus is lost from the Name field.
- The Plate Name field in the Experiment Explorer (“Experiment hierarchy” on page 291) is automatically updated when the new name is entered and validated by the database.
- You can enter up to 50 alpha-numeric characters in the **Name** field.
- Upon validation, the software automatically removes leading and trailing spaces, and converts consecutive spaces to single spaces.
- If you attempt to enter invalid characters, a warning balloon indicates the error condition, and the invalid characters do not appear in the Name field.
- The following characters are not allowed: \ / : \* ? ” < > | %.

- If an invalid name is present in the Name field when focus is lost, the name reverts to the previous valid name.
- Duplicate names are not permitted. If the entered name already exists, a numerical suffix in parentheses is added to the Plate name when the focus is lost; for example, **Plate(2)**. If the addition of the suffix causes the Plate name to exceed the 50-character limit, the entered name is truncated.

## Plate type

- The Plate Type dropdown list allows you to select the type of micro-titer plate (U-, V-, or flat-bottom) that is being used for the current Plate (96- or 384-well). The following options are available:
  - 96 Well Round/U Bottom
  - 96 Well Flat Bottom
  - 96 Well Conical/V Bottom
  - 96 Deep Well Round/U Bottom
  - 96 Deep Well Conical/V Bottom
  - 384 Well Round/U Bottom
  - 384 Well Flat Bottom
  - 384 Well Conical/V Bottom
  - 384 Deep Well Round/U Bottom
  - 384 Deep Well Conical/V Bottom
- You cannot switch between 96- and 384-well plates. The dropdown list is limited to 96- or 384-well plate options based on the original Plate selected when creating the new Experiment as described in “Create a plate experiment” on page 609.

## Options



- **Use Foil Cover** allows you to set up a Plate Experiment using a foil cover to protect the sample plate from condensation or evaporation. When the **Use Foil Cover** option is enabled, the autosampler disables the probe collision sensor, which allows the use of a foil cover on the plate.
- **Use Cooling Block** allows you to set up a Plate Experiment using a cooling block. When selected, the autosampler accounts for the extra height that the cooling block adds to the plate specification.
- **Use Foil Cover** and **Use Cooling Block** options are available only when a CytKick™ Max™ Autosampler is connected to the Attune™ instrument. Otherwise, they are not visible.
- When available, you can also set the **Use Foil Cover** and **Use Cooling Block** options in the New Experiment dialog (“New Experiment dialog” on page 606).

## Plate ID

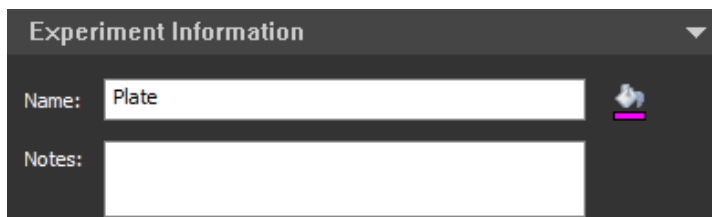
- The Plate ID field allows 50 characters to be entered to identify the selected Plate. Any printable character can be entered.
- The Plate ID is applied to the Plate and is used by any Tubes added to the Experiment.

## Notes

- You can enter up to 500 characters in the Notes field to provide relevant information about the Plate; any printable character is permitted.
- The notes are saved with all Samples in the Plate, including any Tube Experiments added to the Plate.
- Two lines of space are given by default. If you enter more text than can fit in the available space, a vertical scroll bar and the “...” button is displayed adjacent to the text field.
- Clicking the ... **button** opens the *Notes dialog*, which contains an editable text field that allows up to 500 characters to be displayed and edited. You can enter any printable character.

## Experiment information

*Experiment information* options are available when the Heat Map view (Chapter 6, “Heat Map View”) is in focus and any Sample is selected.



## Name

- The **Name** field displays the name of the currently selected Experiment.
- You can modify the name of the selected Experiment by entering the desired name in the **Name** field. The entered name is automatically saved and applied to the Experiment once the focus is lost from the Name field.
- The Experiment Name field in the Experiment Explorer (“Files view” on page 287) is updated automatically when the new name is entered and validated by the database.
- You can enter up to 50 alpha-numeric characters in the **Name** field. Upon validation, the software automatically removes leading and trailing spaces, and converts consecutive spaces to single spaces.
- If you attempt to enter invalid characters, a warning balloon indicates the error condition, and the invalid characters do not appear in the Name field. The following characters are not allowed: \ / : \* ? ” < > | %.

- If an invalid name is present when focus is lost, then the name reverts to the previous valid name.
- Duplicate names are not permitted. If the entered name already exists, a numerical suffix in parentheses is added to the Experiment name when the focus is lost; for example, **Experiment(2)**. If the addition of the suffix causes the name to exceed the 50-character limit, the entered name is truncated.

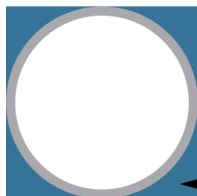
### Experiment color picker

- Clicking the **color icon** opens a standard color picker dialog, which allows you to customize the color for the current Experiment.



- The Experiment color appears on the Heat Map view.

#### Well sample



#### Tube sample



Experiment  
color

### Notes

- The Notes field for the Experiment information works the same way as the Notes field for the Plate information (“Notes” on page 470).

## Group information

*Group information* options are available when the Heat Map view is in focus and any Samples belonging to the same Group are selected.



## Name

- The **Name** field displays the name of the currently selected Group.
- You can modify the name of the selected Group by entering the desired name in the **Name** field. The entered name is automatically saved and applied to the selected Group once the focus is lost from the Name field.
- The Group Name field in the Experiment Explorer (“Files view” on page 287) is updated automatically when the new name is entered and validated by the database.
- You can enter up to 50 alpha-numeric characters in the **Name** field. Upon validation, the software automatically removes leading and trailing spaces, and converts consecutive spaces to single spaces.
- If you attempt to enter invalid characters, a warning balloon indicates the error condition, and the invalid characters do not appear in the Name field. The following characters are not allowed: \ / : \* ? ” < > | %.
- If an invalid name is present when focus is lost, then the name reverts to the previous valid name.
- Duplicate names are not permitted. If the entered name already exists, a numerical suffix in parentheses is added to the Group name when the focus is lost; for example, **Group(2)**. If the addition of the suffix causes the name to exceed the 50-character limit, the entered name is truncated.

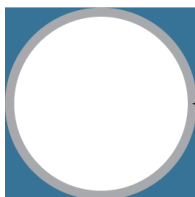
## Group color picker

- Clicking the **color icon** opens a standard color picker dialog, which allows you to customize the color for the current Group.

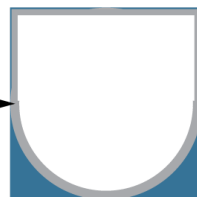


- The Group color appears on the Heat Map view.

### Well sample



### Tube sample



Group  
color

- This option is disabled when the selected Well or Tube is a Compensation control. When multiple Wells are selected, the Group color is not applied to Compensation controls.

## Notes

- The Notes field for the Group information works the same way as the Notes field for the Plate information (“Notes” on page 470).

## Sample information

*Sample information* options are available when the Heat Map view is in focus and a Tube or a Well sample is selected.



- The Name field is hidden if multiple Samples are selected in the Heat Map view.
- The Notes section is enabled and visible when multiple Tube or Well samples are selected.

## Name

- The **Name** field displays the name of the currently selected Sample.
- You can modify the name of the selected Sample by entering the desired name in the **Name** field. The entered name is automatically saved and applied to the Sample once the focus is lost from the Name field.
- The Sample Name field in the Experiment Explorer (“Files view” on page 287) is updated automatically when the new name is entered and validated by the database.
- You can enter up to 50 alpha-numeric characters in the **Name** field. Upon validation, the software automatically removes leading and trailing spaces, and converts consecutive spaces to single spaces.
- If you attempt to enter invalid characters, a warning balloon indicates the error condition, and the invalid characters do not appear in the Name field. The following characters are not allowed: \ / : \* ? ” < > | %.
- If an invalid name is present when focus is lost, then the name reverts to the previous valid name.
- Duplicate names are not permitted. If the entered name already exists, a numerical suffix in parentheses is added to the Sample name when the focus is lost; for example, **Sample(2)**. If the addition of the suffix causes the name to exceed the 50-character limit, the entered name is truncated.

## Notes

- The Notes field for the Sample information works the same way as the Notes field for the Plate information (“Notes” on page 470).

# Customize Overlay options

## Overview

The **Customize Overlay** panel is used for customizing the appearance of **Overlay** and **Gallery plots** in an **Overlay**. The options displayed depend on whether **Overlay plots** or **Gallery plots** are selected.

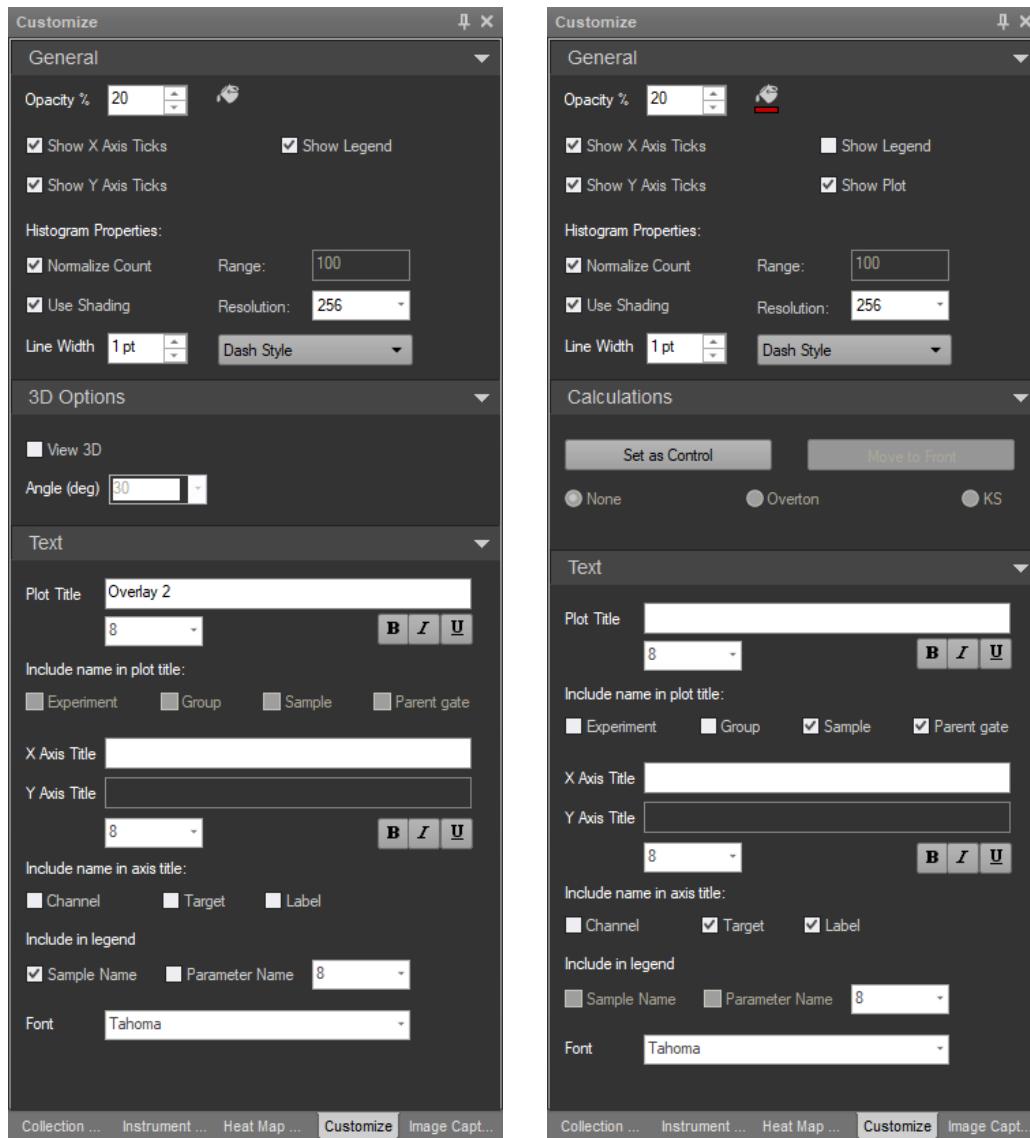


Figure 97 Customize panel when an Overlay plot (left) or a Gallery plot (right) is selected.

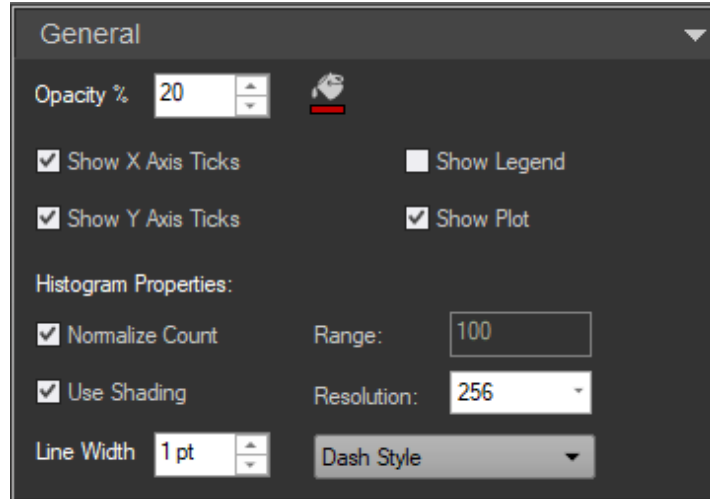
The Customize options are arranged into these functional groups:

- **General** (“General options” on page 475)
- **3D options** (available only for Overlay plots) (“3D Options” on page 479)
- **Calculations** (available only for Gallery plots) (“Calculations” on page 481)
- **Text** (“Text” on page 483)



## General options

**General** options control line style and plot display options. They are available for both **Overlay plots** and **Gallery plots**.



### Opacity

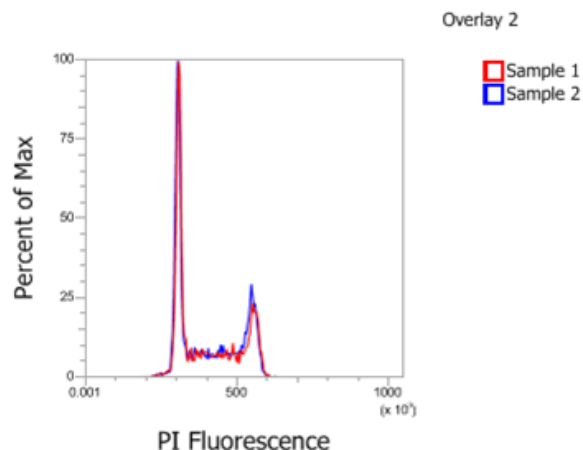
- The **Opacity** option controls the opacity of the line color for Gallery or Overlay Histograms or the data color for dual-parameter plots.
- To change the **Opacity**, enter a number between **20** and **100** into the **Opacity text field**.
- If Histograms are filled, the opacity option applies to both the histogram line and the fill of the histogram.

### Data color

- Click **Data Color** to open the color picker dialog, which allows you to choose the line color for the selected Histograms or the data color for selected dual-parameter plots.
- This option is available when any individual or a combination of gallery plots is selected.
- The selected color is applied to all selected plots.

### Show legend

- Select **Show Legend** to display the legend for an Overlay plot or to display the plot statistics as a legend for a Gallery plot.
- The legend appears to the right of the plot.



### Show X-Axis and show Y-Axis ticks

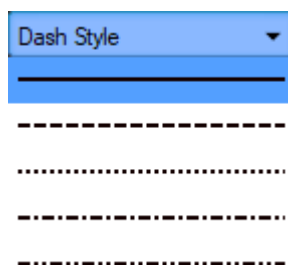
- By default, **Show X-Axis Ticks** and **Show Y-Axis Ticks** options are checked.
- When the **Show X-Axis Ticks** option is checked, the X-axis tick marks are displayed on selected plots.
- When the **Show Y-Axis Ticks** option is checked, the X-axis tick marks are displayed on selected plots.
- When the options are unchecked, the corresponding tick marks are not shown on the selected plots.

### Line width

- The **Line Width** option is only active when Histogram plots in the Gallery or Histogram-based Overlay plots are selected.  
This option is unavailable when only dual-parameter plots are selected.  
If a mixture of plot types are selected (Histograms and dual-parameter plots), this option is available, but only applies to Histogram plots.
- The **Line Width text box** allows adjustment of the line width of selected histogram plots. You can directly enter an integer between 1 and 5 in the textbox.  
Alternatively, you can use the **up** and **down arrows adjacent to the text box** to increase or decrease the line width between **1** and **5**.
- If you enter a line width value that is outside the allowed range, the number is automatically adjusted to the nearest limit.
- If you select different line widths for individual gallery plots, the line width textbox initially appears blank with the edit box active.
- Click the **up arrow** to increase the line width in steps of 1 for each gallery plot and the overlay plot.
- Click the **down arrow** to decrease the line width in steps of 1 for each gallery plot and the overlay plot.
- The up and down arrow buttons have no effect if the edit field is blank.

## Dash Style

- The **Dash Style** option is only available when **Histogram Gallery plots** or **Histogram Overlay plots** are selected.
- The option is not visible when only dual-parameter plots are selected.
- If you select a mixture of plot types (Histograms and dual-parameter plots), this option is visible, but only applies to the selected histogram plots.
- The **Dash Style** dropdown displays a list of available dash styles (solid, dash, dots, dash dot, dash dot dot)
- The dash style selected from the dropdown list applies to all selected Gallery plots or individual members of Overlay plots.
- The default dash type is a solid line.



## Use Shading

- **Use Shading** controls the shading for the selected **Histogram** plots.
- This option applies only to **Histogram Overlay plots** or **Gallery plots**.

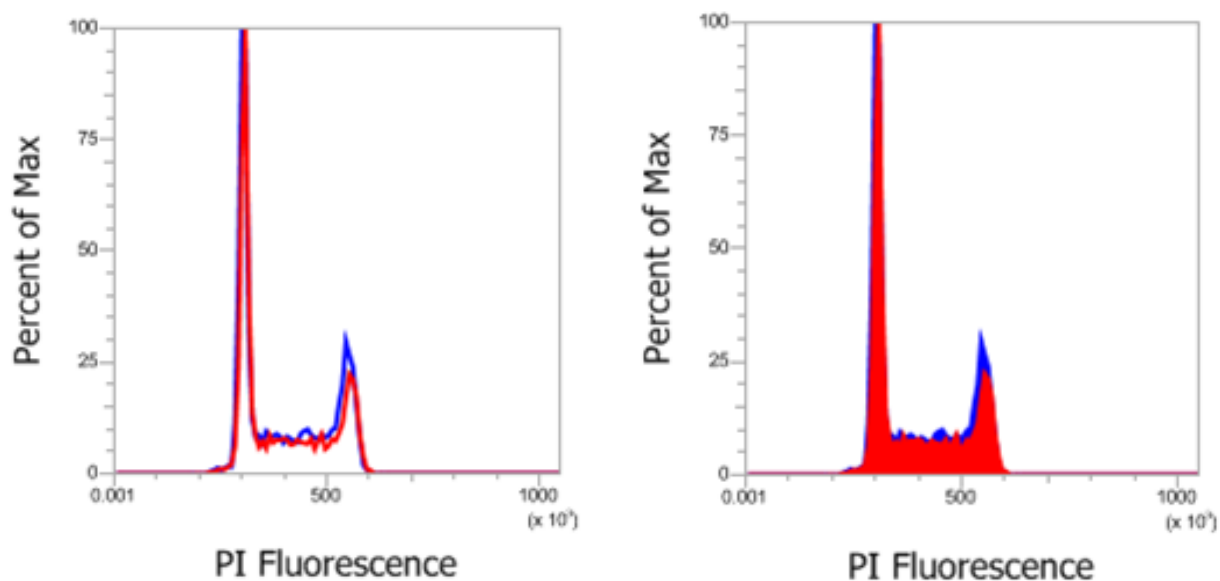


Figure 98 Use Shading - OFF (left) and ON (right)

## Normalize Count

- **Normalize Count** controls the scaling for the selected **Histogram plots**.
- This option applies only to Histogram plots, and it is only displayed when one or more Histogram plots are selected.
- When **Normalize Count** is deselected, the scale label is **Count**.
- When **Normalize Count** is selected, the Y-axis scaling is changed such that the plot is scaled to the Histogram peak value.

The scale displayed is a percentage scale from 0% to 100%, where 100% is the Histogram peak value.

The scale label is **Percent of Max**.

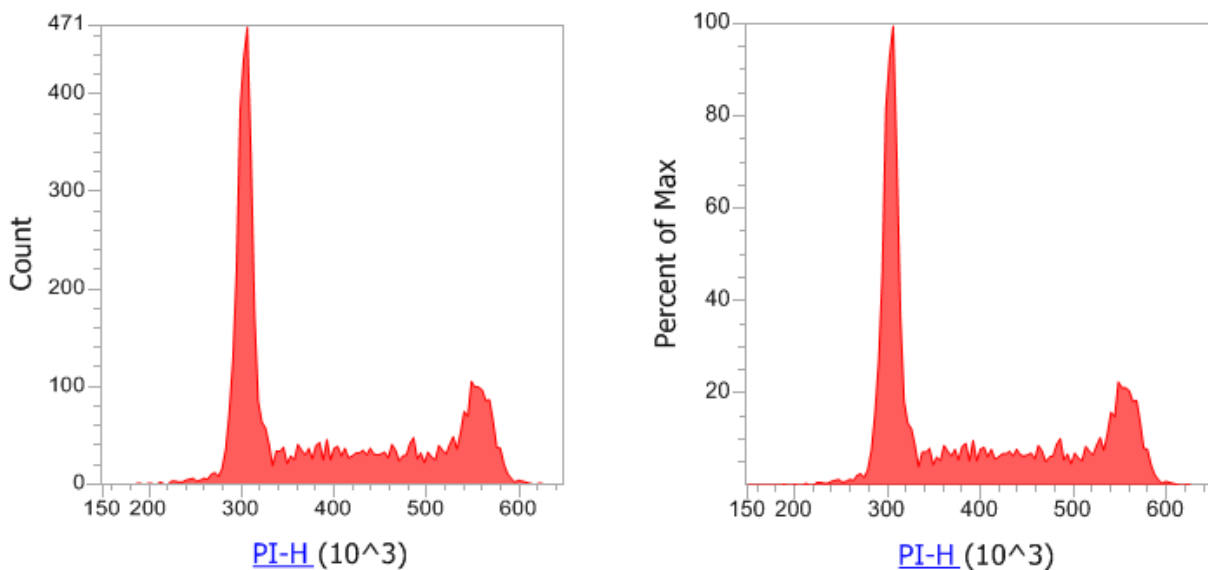


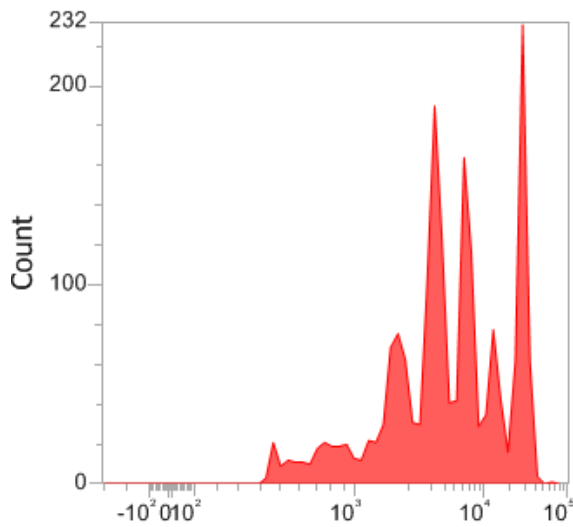
Figure 99 Normalize Count - OFF (left) and ON (right)

## Range

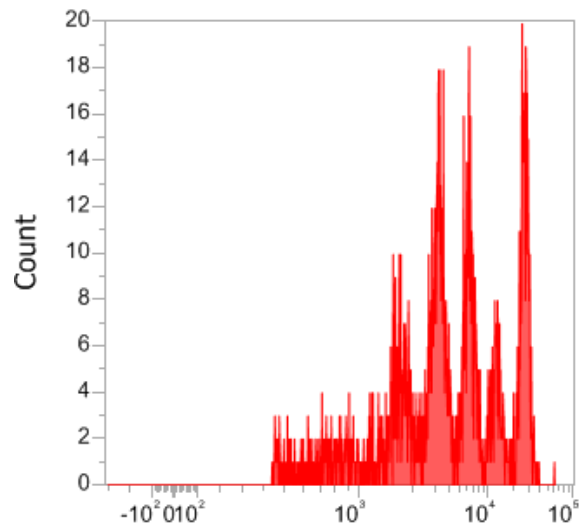
- **Range** sets the upper y-axis range for the selected histograms. The lower end of the range is always set to 0.
- To change the **Range**, enter the desired value in the **Range text field**. Range will not be automatically set and must be manually adjusted.
- The **Range** option is only enabled when the **Normalize Histogram** is deselected.

## Resolution

- **Resolution** determines the resolution for the selected histograms and sets the bin size for drawing the histogram data.
- The available resolutions are **64, 128, 256, 512, and 1024**.
- The default resolution is **1024**.



[8.3 Tg T cells - CellTrace™ Violet-A](#)

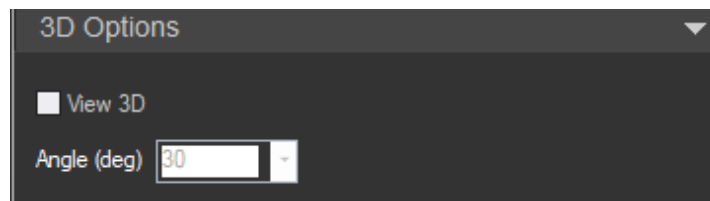


[8.3 Tg T cells - CellTrace™ Violet-A](#)

Figure 100 Resolution: 64 (left) and 1024 (right)

## 3D Options

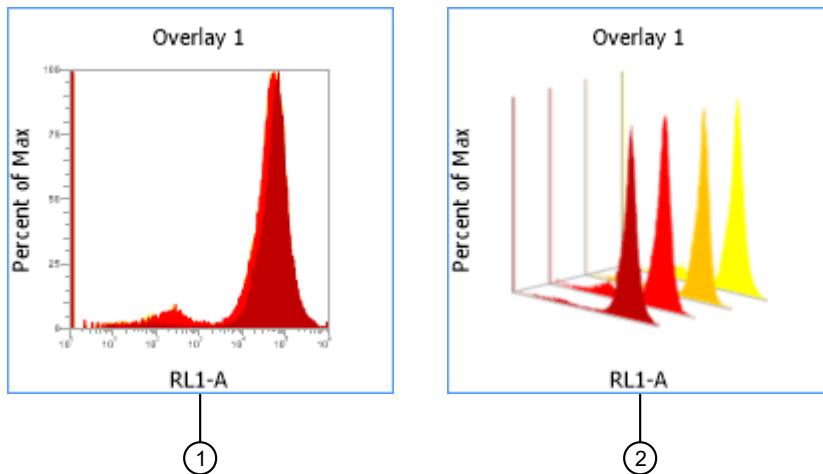
**3D Options** enable you to display **Overlay plots** three dimensionally and set the view option for **3D plots**.



**3D Options** are available only when one or more **Overlay plots** are selected. You cannot view **Gallery plots** in the **3D mode**.

## View 3D

- Select **View 3D** to show selected **Overlay plots** in the **3D mode**.



① Overlay plot in Overlay mode

② Overlay plot in 3D mode

## Angle (deg)

- **Angle (deg)** determines the viewing angle of the **Overlay plot** in the **3D mode**.
- The **Angle (deg)** dropdown becomes active only when **View 3D** is selected.
- **Angle (deg)** dropdown displays a list of angles (**5, 15, 30, 45, 60, 75** degrees). Increasing the angle rotates the 3D plot clockwise.

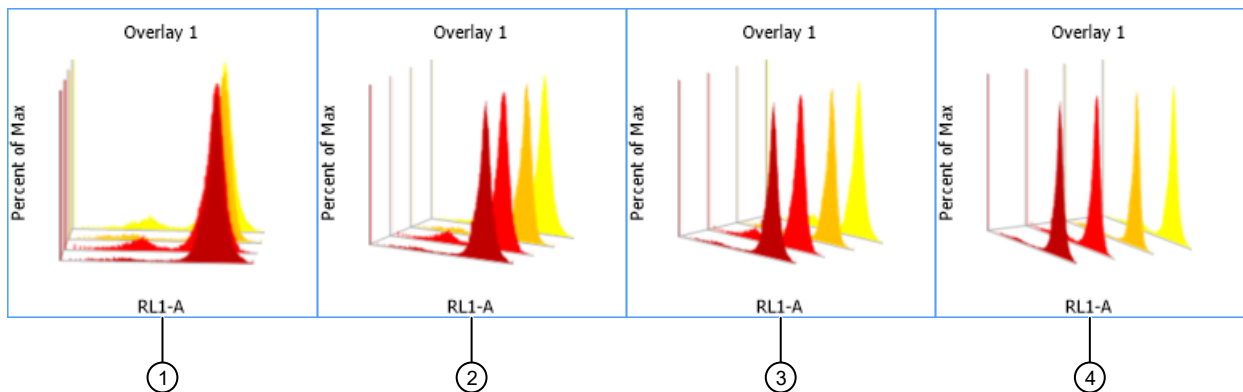


Figure 101 Example 3D plots with viewing angles of 5, 30, 45, and 60 degrees.

① Angle at 5 degrees

② Angle at 30 degrees

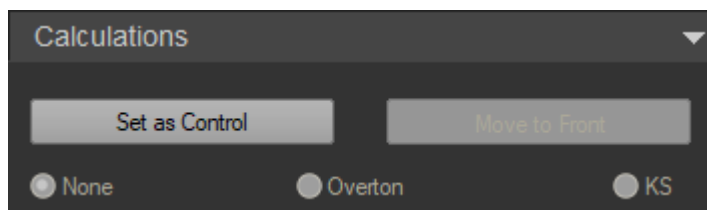
③ Angle at 45 degrees

④ Angle at 60 degrees

## Calculations

The **Calculations** functional group is displayed only for **Gallery plots**.

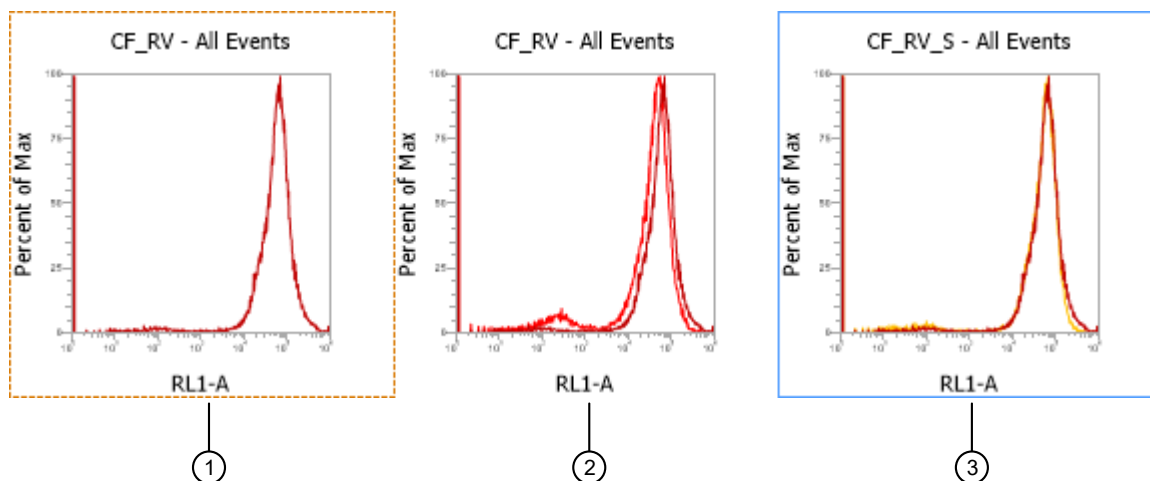
The **Calculations** options are available only for single-parameter plots; they are not displayed if a dual-parameter **Gallery plot** is selected.



### Set as Control

- **Set as Control** option designates the selected plot as the **Control plot** and overlays all **Gallery plots** with it to enable direct visual comparison of several **Test plots** to one **Control plot**.

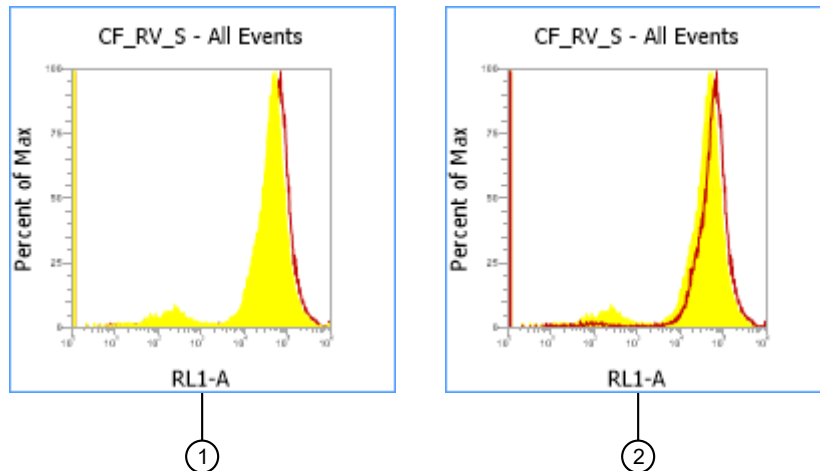
The **Set as Control** button is available only when a single **Gallery plot** is selected. By default, **Set as Control** is deselected.



- ① Control plot
  - ② Gallery plot (Test plot)
  - ③ Gallery plot (Test plot, selected)
- The plot that is set as a **Control plot** is always moved to the first position in the **Gallery** and highlighted in a dashed orange rectangle.
  - To remove a Control plot, click **Remove Control**.

## Move to Front

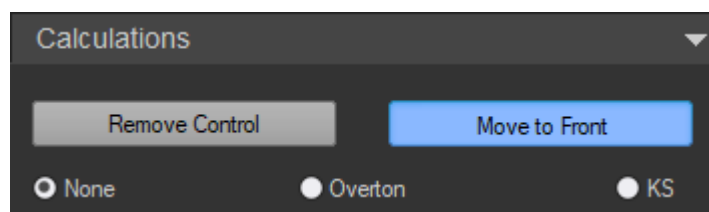
- Click **Move to Front** to show the **Control plot** overlaid on each **Gallery plot** in front of the **Test plot**.  
Click the button again to move the control plot behind the selected plot in the **Overlay gallery**.
- **Move to Front** button is available only when the **Set as Control** option has been selected.
- By default, the **Move to Front** option is not selected and the **Test plot** is displayed in front of the **Control plot**.



- ① Move to Front – OFF: Control plot (red) is behind the Test plot (yellow)  
 ② Move to Front – ON: Control plot (red) is in front of the Test plot (yellow)

## Calculation method options

- **Calculation method** options are only available after a **Gallery plot** has been set as the **Control plot**.



- By default, **None** is selected and no calculations are applied.
- The **Overton** option compares the **Control plot** to each **Gallery plot** using Overton's cumulative statistics calculation ("Overton's cumulative statistics" on page 937).
- The **KS** option compares the **Control plot** and the **Test plot** using the Kolmogorov-Smirnov test ("Kolmogorov-Smirnov test" on page 937) to determine the probability that the two plots are the same.



## Text

**Text** options enable you to customize the plot text on **Overlay** or **Gallery** plots.

The screenshot shows a 'Text' panel with the following controls:

- Plot Title:** A text field containing 'Overlay 1', a font size dropdown set to '8', and buttons for Bold (B), Italic (I), and Underline (U).
- Include name in plot title:** Four checkboxes for 'Experiment', 'Group', 'Sample', and 'Parent gate', all of which are currently unchecked.
- X Axis Title:** A text field.
- Y Axis Title:** A text field, a font size dropdown set to '8', and buttons for Bold (B), Italic (I), and Underline (U).
- Include name in axis title:** Three checkboxes for 'Channel', 'Target', and 'Label', all of which are currently unchecked.
- Include in legend:** Two checkboxes, 'Sample Name' (checked) and 'Parameter Name' (unchecked), followed by a font size dropdown set to '8'.
- Font:** A dropdown menu currently set to 'Tahoma'.

**Text** options consist of the **Text group** and the **Legend group** controls.

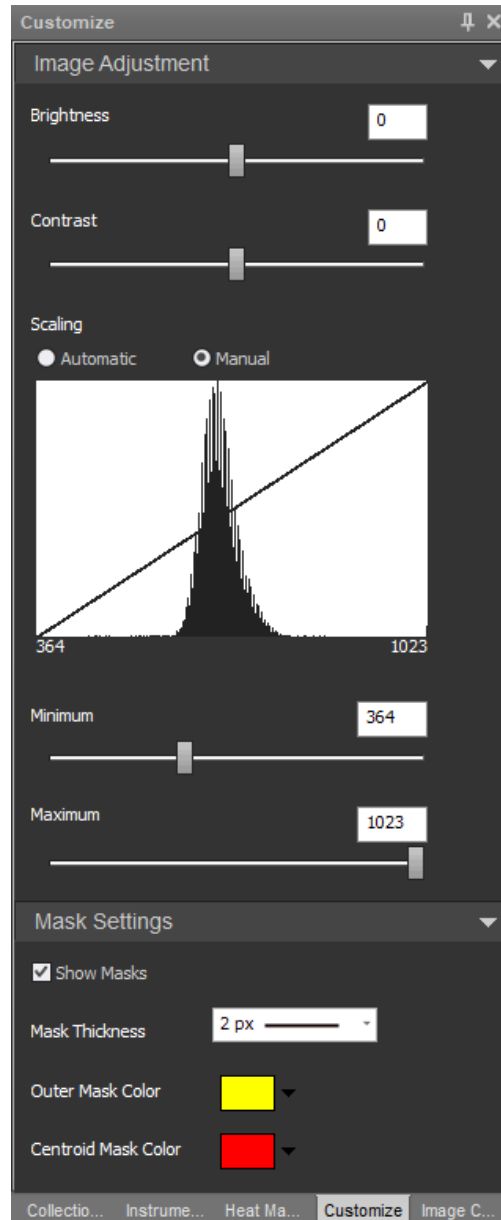
- The **Text group** options are the same as those described for the **Text group** in “Text group (plots only)” on page 442.
- The **Legend group** options contain the **Sample name** and **Parameter name** checkboxes.
- Select the **Sample name** and/or the **Parameter name** checkbox to show the corresponding legend on the plot. You can select either one or both options for display.
- You can edit the legend font size using the **Font size** control. Font size is selected from the **Font size dropdown**, or by entering a number between **6** and **72** into the **font size text field**.

# Customize Image options

## Overview

The **Customize** panel contains the **Image Adjustment** and **Mask Settings** controls, which apply to the active image in the **Image View** tab or in a **Cell Image Container** in the Workspace.

**Note:** The **Customize Image** options are not available for **Workspace images**, which are distinct from images in **Cell Image Containers** or **Image View tab** images.



## Image Adjustment

**Customize** panel **Image Adjustment** group enables you to adjust the **Brightness** and **Contrast** of the cell image in a **Cell Image Container** in the **Workspace** or the active image in the **Image View** tab.

The **Image Adjustment** options are not available for **Workspace images**, which are distinct from images in **Cell Image Containers** or **Image View tab** images.

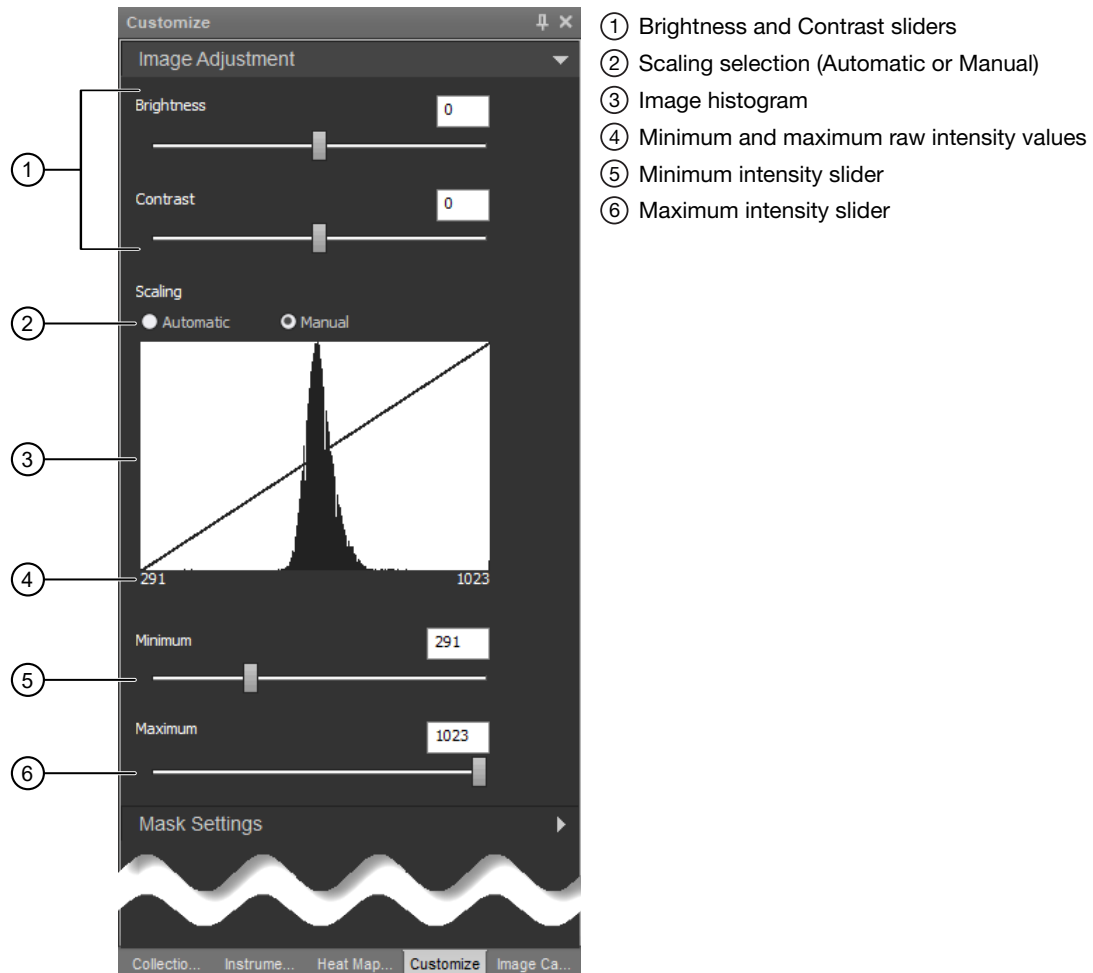


Figure 102 Customize panel Image Adjustment controls

## Brightness and Contrast

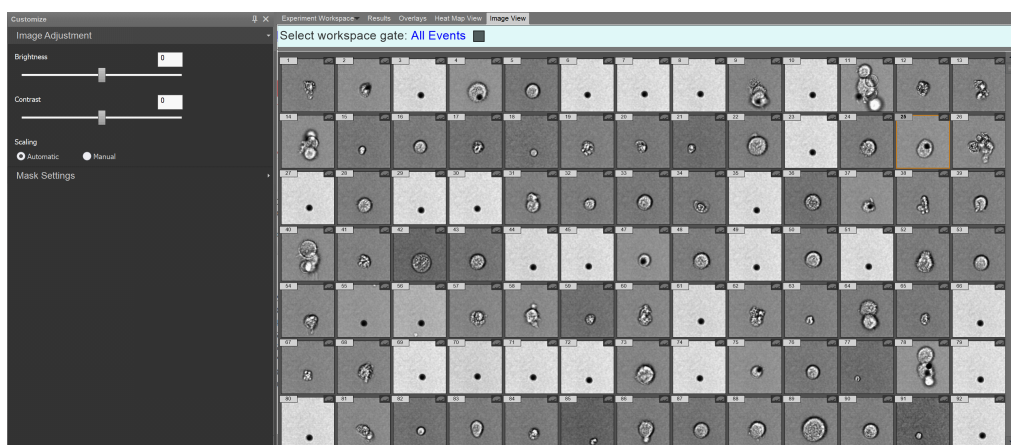
- Slider controls and corresponding text edit fields enable the change of **Brightness** or **Contrast** values.
- Changes to brightness and contrast are applied to all images in the image gallery and view.
- The **Brightness** control allows a value from **-255** to **255**.  
The **Contrast** control allows a value from **-100** to **100**.
- For images in **Cell Image Containers** in the **Workspace**, the **Contrast** and **Brightness** settings are applied on a per cell image basis and the setting persists with the **Workspace**.  
If multiple cell image containers are selected in the **Workspace**, the values displayed in the **Image Adjustment** panel depend on whether the values for all selected images are the same.  
If the brightness and contrast values are different for the selected images, the slider shows the value set at **0** and the edit control is blank for the corresponding setting.  
If the brightness and contrast values match for the selected images, they show the actual value.  
If the value for contrast or brightness is changed, the new value is applied to all selected cell images.
- For images in the active image area and image gallery of the **Image View** tab, the change to brightness and contrast is set globally for all cell images. The brightness and contrast values persist with the **Workspace**.

## Image scaling

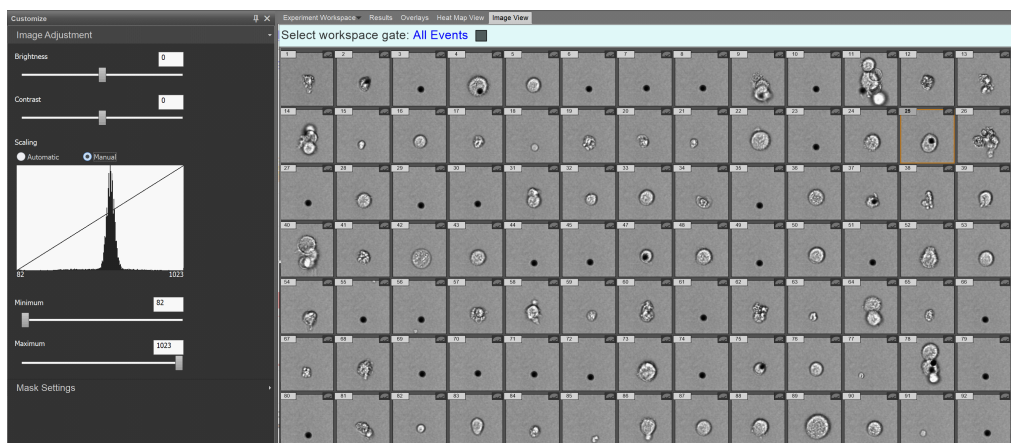
The Attune™ CytPix™ Flow Cytometer captures images with 10 bits of dynamic range (0 to 1023) and stores them as 16-bit greyscale TIFF image files (maximum intensity of 65536). Because computer monitors can only display RGB between 0 to 255 per channel, the range that a single 8-bit byte can offer (which enables 16,777,216 colors –  $256^3$  colors and another 256 levels of opacity), images greater than 8 bits have to be scaled down to 8 bits per channel.

By default, the Attune™ Cytometric Software scales down the images such that the minimum intensity in each image becomes 0 and the maximum intensity becomes 255. This automatic scaling can result in visual artifacts, such as black images or images in which the background appears differently in each image.

The Attune™ Cytometric Software enables you to manually scale the captured images, so that all images have the same minimum and maximum value (that is, they are normalized to a fixed intensity range). Setting a fixed-range normalization value enables images to be compared, where image scaling is identical.



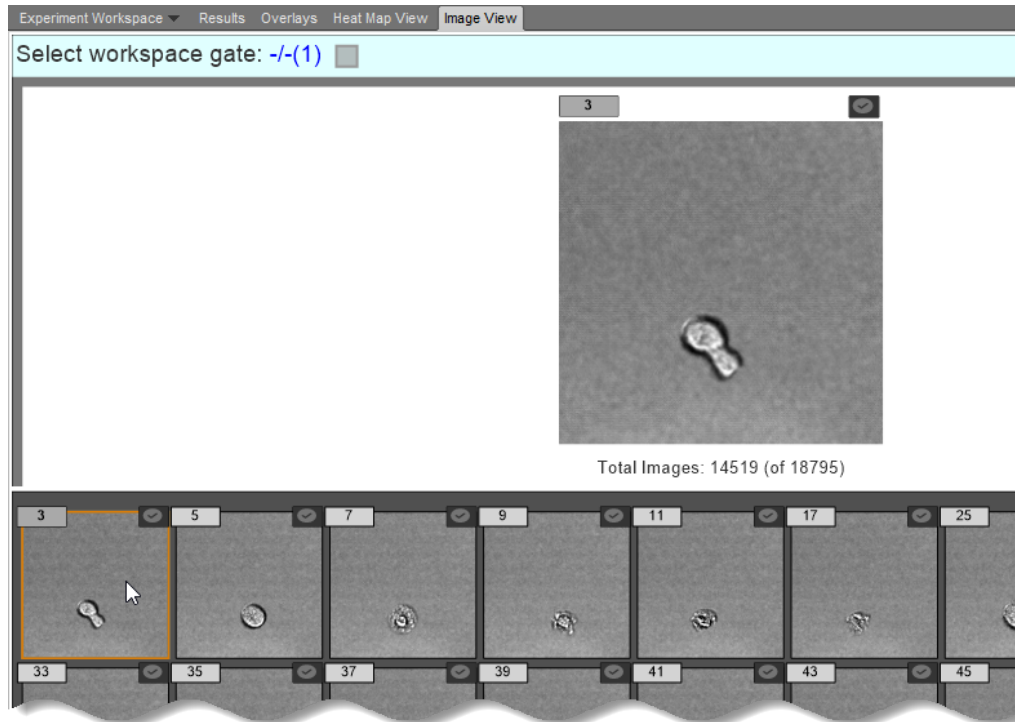
**Figure 103 Automatic scaling** Automatic scaling (default) sets the minimum and maximum intensity within each image to 0 and 255, respectively.



**Figure 104 Manual scaling** Manual scaling enables you to set the minimum and maximum intensity for each image such that the specified minimum and maximum values are scaled to 0 and 255, respectively, which normalizes the images to a fixed intensity range.

## Manually adjust image scaling

1. In **Image View**, select the image to use for manual scaling.



2. In the **Customize** panel, select **Manual** for **Scaling** under **Image Adjustment**.  
The software displays the histogram for the active image that was selected in **Image View**, and the **Minimum** and **Maximum** values reflect the raw intensity values for the selected image.
3. To adjust image scaling, move the **Minimum** and **Maximum** sliders until the you are satisfied with the images displayed in **Image View**. Alternatively, type in the desired **Minimum** and **Maximum** intensity value.

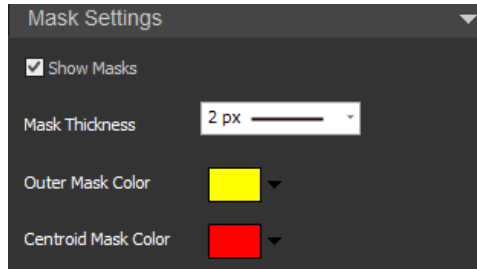
---

**Note:** Image scaling is applied to **all** images in the Image View and gallery, as well as the cell images in the Experiment Workspace. During acquisition, the image scaling is only applied to newly acquired images; the existing preview images remain unchanged until the acquisition completes.

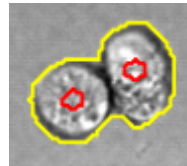
---

## Mask Settings

**Mask Settings** controls enable you to show and hide image masks, and let you to change the line thickness and color of inner and centroid masks.



- When image processing is completed, the software also returns the image masks that were generated by the image processing.



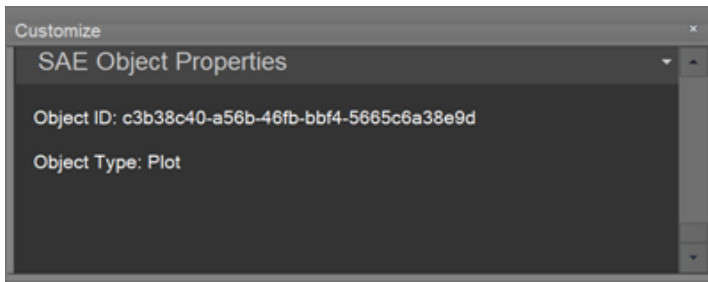
- Image masks are binary representations of an image where an object is identified by the presence of a signal versus its absence. They provide a visual confirmation as to how the image processing “saw” the cell or the particle.
- The pixel coordinates of the mask contour are returned as outputs in the image processing results.
- The results include both the outer masks and the centroid masks (number of spots or cells detected in the Field of View). In the example above, the outer masks are depicted in yellow and the centroid masks in red.
- To show or hide the image masks, select or deselect **Show Masks**.
- To change the line thickness of the mask contour, select the desired thickness in pixels from the **Mask Thickness** dropdown.
- To change the color of the outer or the centroid mask, select the desired color from the **Outer Mask Color** or the **Centroid Mask Color** dropdown.

# SAE object properties

## Overview

When the Attune™ Cytometric Software is in the SAE mode, you can view the ID of workspace objects in the object's customize panel. Each customize panel will have an additional group where the “SAE Object Properties” for the selected workspace object are displayed.

- Each workspace object has a unique ID or GUID. This ID is an immutable property and cannot be changed. This ID is used by the SAE console to enable auditing of workspace objects.
- The workspace objects that have a corresponding GUID are workspace plots, workspace gate, workspace text boxes, workspace statistics boxes, workspace images, overlay plots, and overlay/gallery plots.
- The SAE Object Properties displays the Object ID and Object Type.
- You can copy the Object ID to the clipboard by double clicking on the text of the Object ID.





## Overview

**Image Capture Settings** panel lets you adjust settings that affect the capture of images.

- **Image Capture Settings** panel is only available when using an Attune™ CytPix™ Flow Cytometer.
- By default, the panel is docked to the left of the **Main Application Area**. The docking properties of the panel are described on “Docking locations” on page 58.
- The panel contents are organized into the these groups:
  - **Settings options** (“Settings options” on page 493)
  - **Capture Settings** (“Capture Settings” on page 494)
  - **Camera Settings** (“Camera settings” on page 497)
  - **Image Size and Position Settings** (“Image Size and Position Settings” on page 499)
  - **Advanced** (“Advanced” on page 502)
- Each group can be collapsed and expanded, except **General** panel properties.  
By default, the groups are all expanded, unless otherwise specified below. The expanded or collapsed state persists on a per user basis.
- During Run mode, image capture settings can be adjusted to optimize image capture, and they can be stored at the Sample or Experiment level.
- During recording or after data are recorded, the **Image Capture Settings** are disabled to reflect the state in which the sample was recorded.

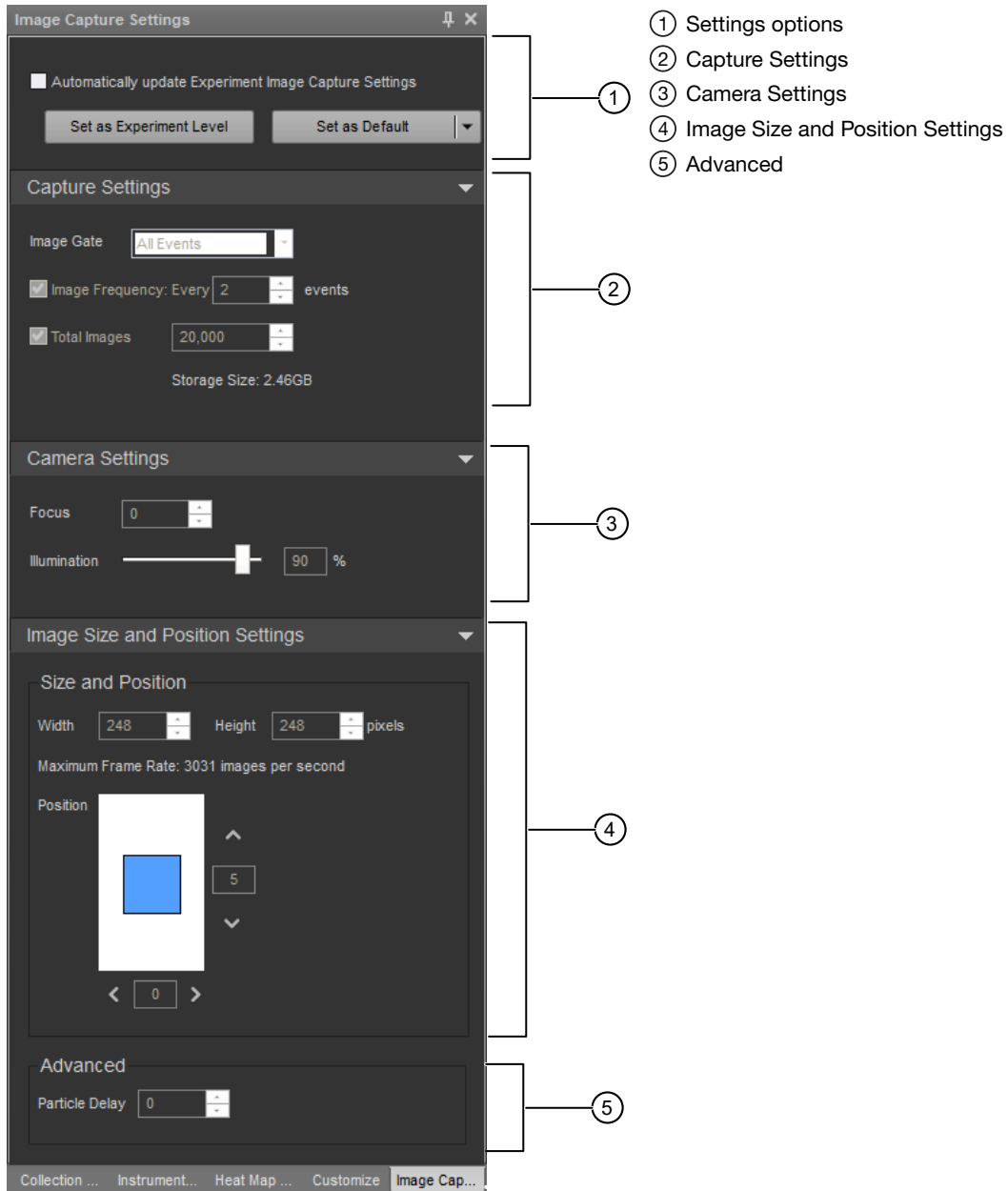
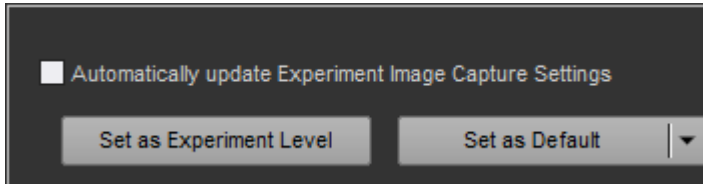


Figure 105 Image Capture Settings panel

## Settings options

**Settings options** section of the **Image Capture Settings** panel contains three controls:



- **Automatically Update Experiment Image Capture Settings** checkbox
- **Set as Experiment Level** button
- **Set As Default** split button

### Automatically update

When selected, **Automatically Update Experiment Image Capture Settings** option automatically applies the changes to **Capture**, **Camera**, and **Image Size** and **Position Settings** to the Experiment-level settings.

This checkbox is disabled if the current active Experiment Explorer item is a **Compensation node** or a **Sample** with **Sample-level Image Capture Settings**.

### Set as Experiment Level

**Set as Experiment Level** updates the appropriate Experiment-level Image Capture Settings with the settings of the active sample (Tube or Well).

Set all applicable Samples (tube or well) in the Experiment to use the new Experiment-level Image Capture Settings.

There is an Experiment-level Image Capture Settings for tubes and another one for plates.

All newly created Samples use the Experiment-level Image Capture Settings until they are activated or edited, at which point a Sample-level Image Capture Setting is created for the applicable samples. This newly created setting is a copy of the Experiment-level Image Capture Settings.

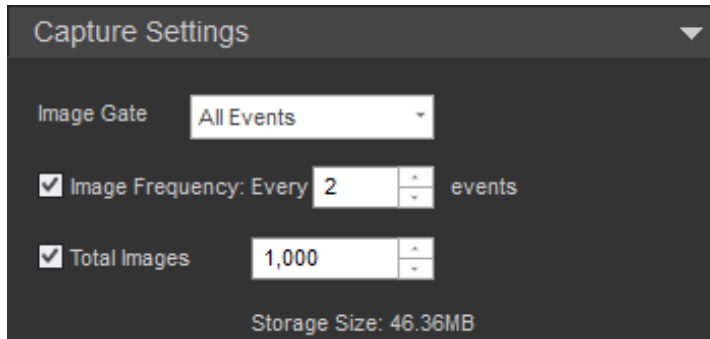
### Set as Default

**Set as Default** split button contains three options:

- **Set as Default** sets the current **Image Capture Settings** as the default for all future Plate or Tube experiments run by the current user.
- **Load** opens the standard **File Open (Import)** dialog, which enables you to import image capture settings from a saved **Image Capture Settings file (\*.aics)** and update the current settings with the imported settings.
- **Export** opens the standard **File Save (Export)** dialog, which enables you to export the current **Image Capture Settings** with a user defined name. If the name selected for the capture settings matches the name of another capture setting in the database, you are prompted for confirmation before the existing image capture setting is overwritten.

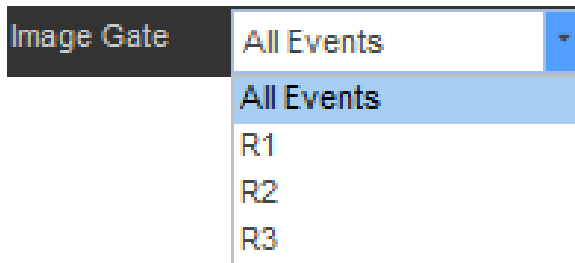
## Capture Settings

**Capture Settings** enables you to set image capture criteria for the sample and contains the **Image Gate**, **Image Frequency**, and **Total Images** controls. It also provides information about the **Storage Size** needed for the captured images based on the capture criteria.



### Image Gate

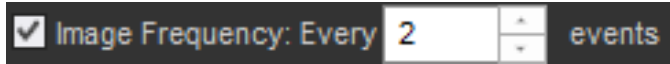
**Image Gate** dropdown enables you to set an image capture gate, so that the images are captured only from a specific subset of events on a **Workspace plot**. Events that are outside of the **Image Gate** are excluded from image collection.



- **Image Gate** dropdown only allows **Rectangle gates** on **dual-parameter plots** and **Histogram gates** on **single-parameter plots** to be used as the **image capture trigger gate**. The default **trigger gate** is **All Events**.
- The coordinates of the **trigger gate** and the corresponding parameter define the image threshold that triggers the system to capture an image if a cell's signal falls within the trigger window's minimum and maximum values (inclusive).
- The image trigger window persists on a per Sample or per Experiment basis.
- If the **trigger gate** is moved during acquisition, the new trigger threshold values are sent to the instrument.  
If the **trigger gate** is deleted during acquisition, the trigger window defaults to **All Events**. In this case, any event triggers image acquisition.
- The image settings persist as part of the Experiment on a per Experiment or Sample basis.
- **Image Gate** does not allow you to set up population hierarchies; it only supports a top-level perspective.

## Image Frequency

**Image Frequency** sets the image capture frequency such that every Nth event that exceeds the imaging trigger setting is imaged. This enables you to capture only a subset of reference images rather than imaging all events and reduces the amount of data storage needed for image collection.



- By default, **Image Frequency** setting is unchecked.
- By default, **Image Frequency** is set to **every 10 events**.
- You can set an **Image Frequency** value of between **every 1** and **10,000 events**.
- If the **Image Frequency** value is changed during a run, the new frequency setting is applied only to new events and does not affect any previously acquired images.
- During data recording or after the recording is completed, the **Image Capture Settings** are disabled to reflect the state in which the Sample was recorded.

## Total Images and Storage Size

**Total Images** sets the maximum number of images to be acquired during a run (if selected). This lets you limit the number of images collected and reduces the amount of data storage needed for image collection.



- By default, **Total Images** is checked and set to **10,000 events**.
- You can set a **Total Images** value of between every **1** and **30,000 events**.
- If the **Total Images** value is changed during a run, the new setting is applied to the runtime total. If the **Total Images** value is changed such that the total number of imaged events is less than the number of currently captured events, no new images are acquired.
- During data recording or after the recording is completed, the **Image Capture Settings** are disabled to reflect the state in which the sample was recorded.

**Storage Size** shows the amount of storage needed to store the captured images when the **Total Images** option is enabled.

- The **Storage Size** value is calculated from the total number of images and the image size (based on width and height specified in **Image Size** and **Position Settings**).
- If the number of images, image width, or image height are changed, the **Storage Size** is automatically recalculated and the new value is displayed.
- When the **Total Images** option is disabled, **Estimated Total Images** is displayed instead of **Storage Size**.

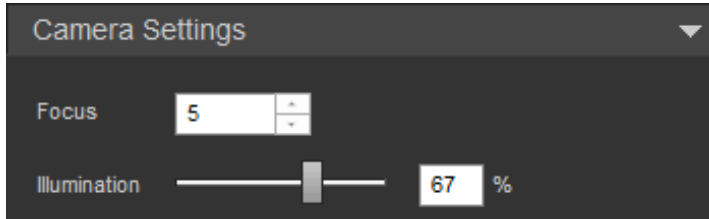
---

**Note:** By default, captured images are stored in the following location:  
C:\Users\Public\Documents\Life Technologies\AttuneCytPix\_Images. Users with Administrator or System Administrator roles can select a different image storage path using the Attune™ Database Utility (Chapter 29, “Attune™ Database Utility”).

---

## Camera settings

**Camera Settings** lets you control the camera **Focus** and **Illumination**.



### Focus

**Focus** enables you to adjust the focus of the camera used for image capture.



- Camera focus is preset to the optimal position with the default focus value of **0**.
- To change the focus of the camera, enter a number into the **Focus edit field** or adjust the value using the adjacent **up** and **down** controls.
- The **up** and **down** controls change the focus value by **1**.
- The **Focus edit field** only accepts integers between **-1000** and **1000**.
- If the focus value that was entered exceeds the limits, the focus edit field reverts to the last good value and a warning message instructs you to enter a number within the acceptable range.
- During recording or after data are recorded, the camera settings are disabled to reflect the state in which the sample was recorded.

---

**Note:** Although the focus is preset to the optimal position by service engineers, depending on cell size, cell type, or flow rate, you might select to modify the **Focus** value to suit the needs of your experiment. Usually, a change of **± 10 units** is sufficient to find the ideal focus position for most applications.

---

## Illumination

**Illumination** slider bar and edit field allow you to adjust the camera illumination used for image capture.



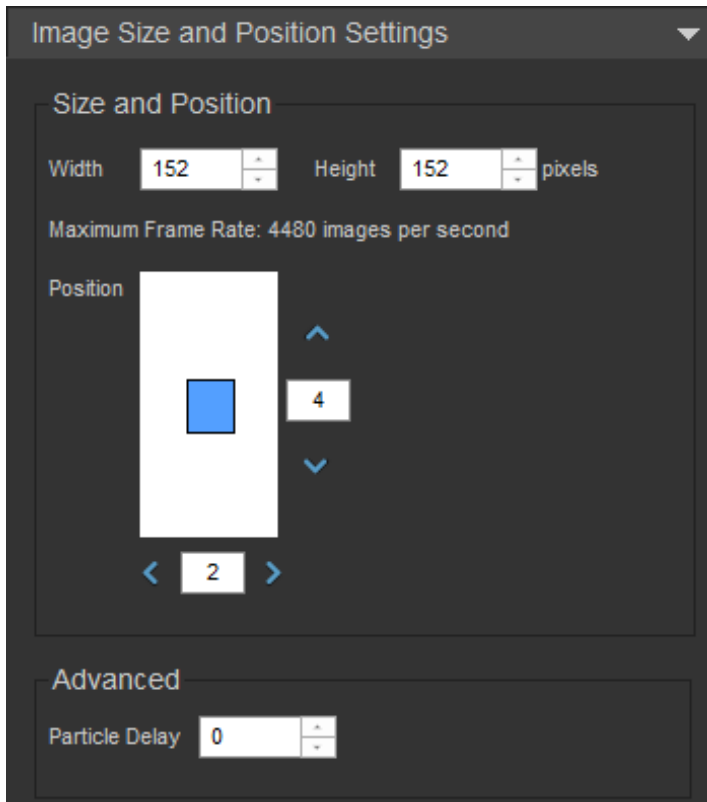
- The default value for camera **Illumination** is **100%**.
- To change the camera **Illumination**, enter a number into the **Illumination edit field** or adjust the value using the **Illumination slider**.
- The **Illumination edit field** only accepts integers between **1** and **100**.  
Clicking the keyboard **up** (or **right**) and **down** (or **left**) **arrow buttons** when the **edit control** is active changes the illumination value by **5 units**.
- The **Illumination slider** has a resolution of **1** unit.  
Clicking the keyboard **up** (or **right**) and **down** (or **left**) **arrow buttons** when the **slider control** is active changes the illumination value by **1 unit**.  
Clicking the keyboard **Page Up** and **Page Down** buttons when the **slider control** is active changes the illumination value by **5 units**.
- If the entered focus value exceeds the limits, the illumination edit field reverts to the last good value and a warning message instructs you to enter a number within the acceptable range.
- During recording or after data are recorded, the **Camera Settings** are disabled to reflect the state in which the sample was recorded.



## Image Size and Position Settings

**Image Size and Position Settings** enable you to adjust the **size** and **position** of the **image window** and to set **particle delay**. It also displays the **image frame rate**, which is calculated based on the height and width of the image window.

During recording or after data are recorded, the **Image Size and Position Settings** are disabled to reflect the state in which the sample was recorded.



---

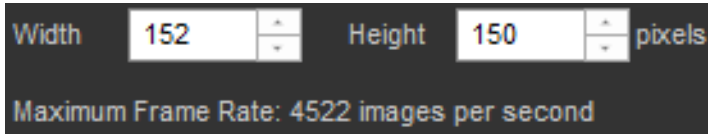
**Note:** Changing image capture settings is not instant; it takes 1–2 seconds for the camera to be cleared and streaming to restart. We do **not** recommend changing the image size and position incrementally with mouse clicks.

---

## Size and Position

### Image Size

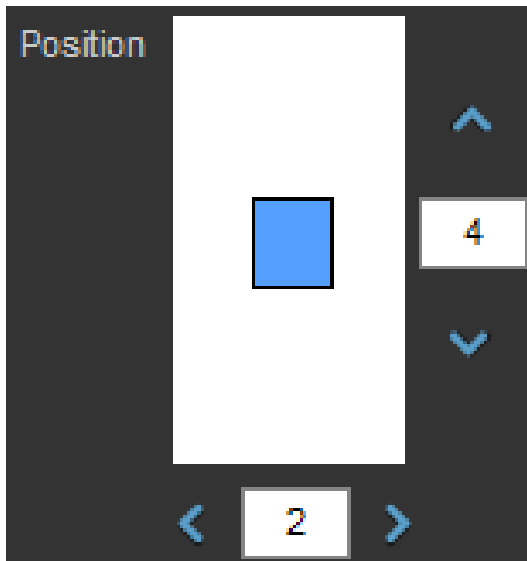
**Image Size** controls consist of **Width** and **Height** controls, which enable you to set the image dimensions.



- **Width** edit field lets you enter a value from **96 pixels** to **248 pixels** for the width of the image window. The **up** and **down** buttons adjust the width by **8 pixels** each click.
- **Height** edit field lets you enter a value from **96 pixels** to **248 pixels** for the height of the image window. The **up** and **down** buttons adjust the height by **2 pixels** each click.
- The default values for **Width** and **Height** are **152 pixels**, which is the optimal size for the image window.
- If you change the image size and capture new images, all the images shown in Image Gallery will have the new aspect ratio as defined by new width and height.
- **Maximum Frame Rate** displays the maximum rate by which the system can trigger the camera for image capture, which is calculated based on the pixel size, image window, and system electronics.

## Position

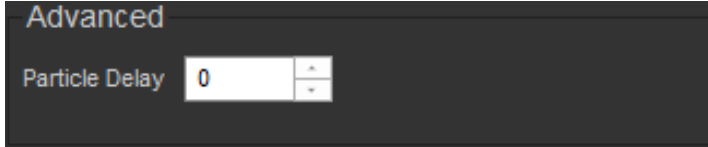
Image **Position** controls allow you to adjust the position the image window within the scanning region. The controls consist of a blue rectangle that represents the **image window** within the **scanning region** and **directional arrow controls** and **position edit fields** to adjust the coordinates of the **image window**.



- The **blue rectangle** represents the **image window** and the **white rectangle** represents the **scanning region** that surrounds it.  
Changing the size of the **image window** using the **Width** and **Height** controls changes the size of the **blue rectangle** that represents it.
- The maximum and minimum size of the **image window** is **248 × 248 pixels** and **96 × 96 pixels** (width × height), respectively.  
96 × 96 pixels correspond to a field of view that is 29 × 29 μm and 248 × 248 pixels correspond to a field of view that is 74 × 74 μm.  
The default size of the **image window** is **152 × 152 pixels** (46 × 46 μm field of view), which is also the optimal size for image collection.
- The size of the **scanning region** is fixed at **450 × 750 pixels** (width × height). The **scanning region** adheres to the coordinate system where the coordinates are (-/-) on the top left and (+/+) on the bottom right.
- By default, the **image window** (blue rectangle) is centered around the coordinates **0** and **0**.
- You can change the position of the **image window** within the **scanning region** using one of the these methods:
  - Drag and drop the **blue rectangle** in a new position within the **scanning region**.
  - Enter the **new position coordinates** into the **X** and **Y coordinate fields**.
  - Use the **directional arrows** to adjust the position of the **image window** by **1 pixel** increments.

## Advanced

**Advanced** group consists of the **Particle Delay** control, which enables you to adjust the image delay relative to the trigger event.



- By default, **Particle Delay** value is set to **0**.
- To change the **Particle Delay**, enter a number into the **Particle Delay edit field** or adjust the value using the adjacent **up** and **down** controls.
- The **up** and **down** controls change the focus value by **1**.
- The **Particle Delay edit field** only accepts integers between **-100** and **100**.

---

**Note:** **Particle Delay** changes the position of the particle to keep it within the **image capture region**. **Size** and **Position** settings move the image capture region within the **scanning area**. If the **Particle Delay** and **Size** and **Position** settings are mismatched, you will not be able to see and image the particle.

To reset the **Position** and **Particle Delay** settings, enter **0/0** for **Position** to center the image field in the scanning region and enter **0** for **Particle Delay**.

---

# 17

## FCS information panel

### Overview

*FCS Information* panel displays the FCS metadata (the text segment portion) of the selected Sample.

**FCS file information**

Filename: 4 COLOR.fcs  
 Filepath: C:\Users\Public\Documents\Life Technologies\AttuneM

---

**Sample Information**

Filename (\$FIL) Sample.fcs  
 Date (\$DATE) 17-Mar-2015  
 Start Time (\$BTIM) 15:31:12  
 End Time (\$ETIM) 15:31:42  
 Institution (\$INST) NA  
 Investigator (\$EXP) NA  
 operator (\$OP) NA  
 Originality (\$ORIGINA... Original  
 Experiment (\$PROJ) concentration  
 Specimen (\$SMNO) LYSED WHOLE BLOOD  
 Sample Source (\$SRC) NA  
 Sample Information (\$... NA  
 Flow Rate (#FLOWRA... 100  
 Comment (\$COM) NA  
 Volume (\$VOL) 50000  
 Timestep (\$TIMESTEP) 0.001  
 Trigger 1 (#TR1) FSC,150000  
 Lost Events (\$LOST) 0  
 Total Events (\$TOT) 16191

---

**Parameters (13)**

Param#	1	2	3	4	5
Name	Time	FSC-A	SSC-A	BL3-A	YL
Stain	Time	FSC-A	SSC-A	cd19-PER...	CD
Target	NA	-	-	cd19	CD
Label	NA	-	-	PERCP-C...	PE
Voltage	NA	370	400	450	310
Range	67108864	1048576	1048576	1048576	104
Bits	32	32	32	32	32
Excitation Wavelength...	NA	NA	NA	NA	NA
Optical Filter	NA	NA	NA	NA	NA
Amplification	0,0	0,0	0,0	0,0	0,0

- The panel is set as a floating panel by default (i.e., its “Dockable” attribute deselected), but it can be set as a dockable panel as described on “Title bar context menu” on page 62.
- The panel header displays the name of the selected Sample, and the complete file path is shown under the panel header.
- The panel contents are organized into the following groups:
  - *Sample information* (“Sample information group” on page 505)
  - *Parameters* (“Parameters group” on page 507)
  - *Compensation* (“Compensation group” on page 508)
  - *System information* (“System information group” on page 509)
  - *Data format* (“Data format group” on page 510)
  - *Other* (“Other group” on page 511)
- Each group can be collapsed and expanded. By default, the groups are all expanded, unless otherwise specified below. The expanded or collapsed state persists on a per user basis.
- The metadata are displayed in a table where the first column is the FCS keyword and the second column is the keyword value.
- Only keywords and keyword values that are present within the FCS file are shown.
- If the panel contents exceed the visible area, vertical and horizontal scroll bars are shown as needed.

---

**Note:** The FCS keyword and keyword values are as described in Appendix A, “FCS file reference”.

---

## FCS file information

### Sample information group

The *Sample information group* contains keywords that describe the FCS file.

The keywords displayed in the Sample information group are listed in the following table and are shown in the sort order indicated in the table.

Sample Information	
Filename (\$FIL)	Sample.fcs
Date (\$DATE)	17-Mar-2015
Start Time (\$BTIM)	15:31:12
End Time (\$ETIM)	15:31:42
Institution (\$INST)	NA
Investigator (\$EXP)	NA
operator (\$OP)	NA
Originality (\$ORIGINA...)	Original
Experiment (\$PROJ)	concentration
Specimen (\$SMNO)	LYSED WHOLE BLOOD
Sample Source (\$SRC)	NA
Sample Information (\$...)	NA
Flow Rate (#FLOWRA...)	100
Comment (\$COM)	NA
Volume (\$VOL)	50000
Timestep (\$Timestep)	0.001
Trigger 1 (#TR1)	FSC,150000
Lost Events (\$LOST)	0
Total Events (\$TOT)	16191

Sort	Keyword	Description
1	\$FIL	Filename
2	GUID – BD™ Keyword	GUID
3	\$DATE	Date
4	\$BTIM	Start Time
5	\$ETIM	End Time
6	\$INST	Institution
7	\$EXP	Investigator

(continued)

Sort	Keyword	Description
8	\$OP	Operator
9	\$LAST_MODIFIER	Last Modified By
10	\$LAST_MODIFIED	Last Modified
11	\$ORIGINALITY	Originality
12	\$PLATENAME	Plate Name
13	\$PLATEID	Plate ID
14	\$WELLID	Well Location
15	\$PROJ	Experiment
16	\$SMNO	Specimen
17	\$SRC	Sample Source
18	\$CELLS	Sample Information:
18	#FLOWRATE	Flow Rate
19	\$COM	Comment
19	\$VOL	Sample Volume
20	#TOTALVOLUME	Sample Volume
21	\$TIMESTEP	Timestep
22	\$TR	Trigger
23	\$ABRT	Aborted Coincident Events
24	\$LOST	Lost Events
25	\$TOT	Total Events



## Parameters group

The *Parameters group* contains the parameter description keywords (e.g., \$PnN, \$PnS, etc.) for each parameter.

The number of parameters (Parameters (\$PAR)) is indicated above the table displaying the Parameter keywords.

Param#	1	2	3	4	5
Name	Time	FSC-A	SSC-A	BL3-A	YL
Stain	Time	FSC-A	SSC-A	cd19-PER...	CD3
Target	NA	-	-	cd19	CD3
Label	NA	-	-	PERCP-C...	PE
Voltage	NA	370	400	450	310
Range	67108864	1048576	1048576	1048576	104
Bits	32	32	32	32	32
Excitation Wavelength...	NA	NA	NA	NA	NA
Optical Filter	NA	NA	NA	NA	NA
Amplification	0,0	0,0	0,0	0,0	0,0

The column headers indicate the parameter number and the row labels correspond to the Parameter description keywords as described in the following table.

Sort	Keyword	Description
1	\$PnN	Channel Name
2	\$PnS	Stain
3	#PnTarget	Target
4	#PnLabel	Label
5	\$PnV	Voltage
6	\$PnR	Range
7	\$PnG	Gain
8	\$PnB	Bits
9	\$PnL	Excitation Wavelength
10	\$PnO	Excitation Power
11	\$PnF	Optical Filter
13	\$PnE	Amplification
14	\$PnP	% Emitted Light
14	\$PnT	Detector Type

(continued)

Sort	Keyword	Description
15	\$PnD	Visualization Scale
16	\$PnCALIBRATION	Calibration

## Compensation group

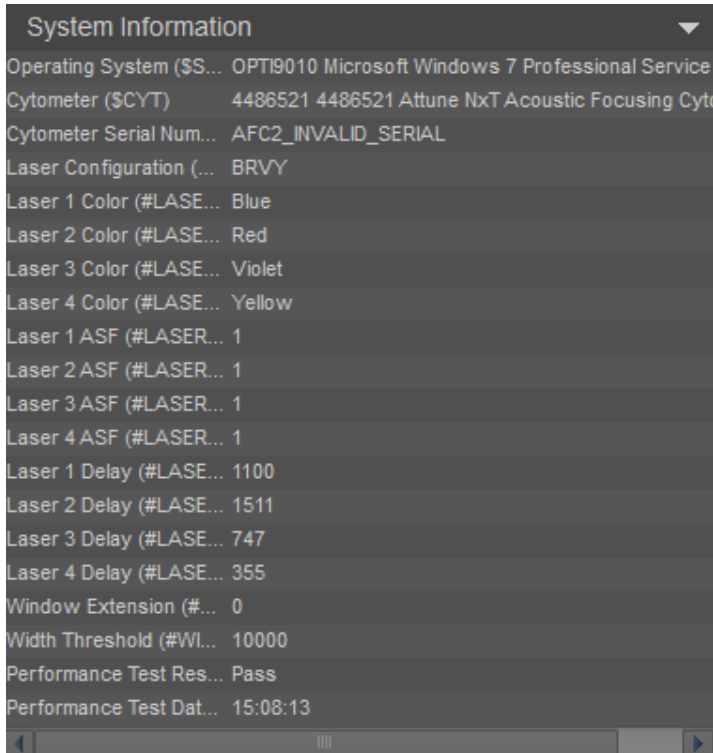
The *Compensation group* displays the spillover values contained within the FCS file in a matrix view where the row and column headers are named using the \$PnN values (i.e., Channel name).

The keywords used to establish compensation (\$Spillover, \$Comp, Spill, Spillover, None) are shown next to the compensation group name.

	BL3-A	YL1-A	RL1-A	VL2-A	BL3-H	YL1-H	RL1-H	VL2-H
BL3-A	100.00	0.04	3.20	0.02	0.00	0.00	0.00	0.00
YL1-A	63.74	100.00	0.07	0.04	0.00	0.00	0.00	0.00
RL1-A	2.48	0.01	100.00	0.01	0.00	0.00	0.00	0.00
VL2-A	0.30	0.01	0.08	100.00	0.00	0.00	0.00	0.00
BL3-H	0.00	0.00	0.00	0.00	100.00	0.00	0.00	0.00
YL1-H	0.00	0.00	0.00	0.00	63.74	100.00	0.00	0.00
RL1-H	0.00	0.00	0.00	0.00	2.48	0.00	100.00	0.00
VL2-H	0.00	0.00	0.00	0.00	0.30	0.00	0.00	100.00

## System information group

The *System information group* contains keywords that pertain to the system. For all custom keywords, keywords without the “#” sign will also be recognized.



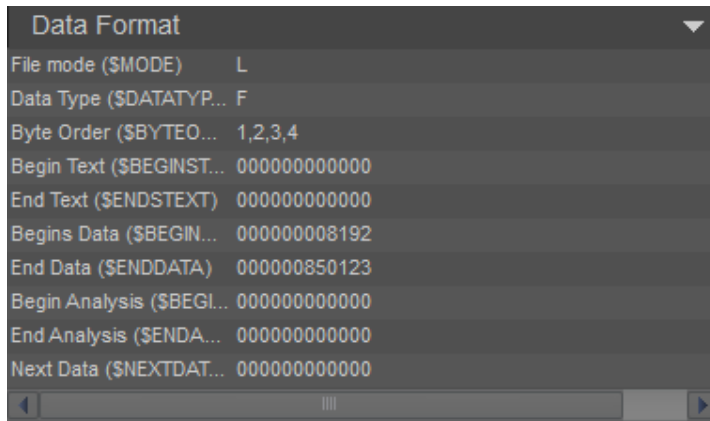
Sort	Keyword	Description
1	\$SYS	Operating System
2	\$CYT	Cytometer
3	\$CYTSN	Cytometer Serial Number
4	#LASERCONFIG	Laser Configuration
5	#LASER1COLOR	Laser 1 Color
6	#LASER2COLOR	Laser 2 Color
7	#LASER3COLOR	Laser 3 Color
8	#LASER4COLOR	Laser 4 Color
9	#LASER1ASF	Laser 1 ASF
10	#LASER2ASF	Laser 2 ASF
15	#LASER3ASF	Laser 3 ASF
16	#LASER4ASF	Laser 4 ASF

(continued)

Sort	Keyword	Description
17	#LASER1DELAY	Laser 1 Delay
18	#LASER2DELAY	Laser 2 Delay
18	#LASER3DELAY	Laser 3 Delay
19	#LASER4DELAY	Laser 4 Delay
19	#WINEXT	Window Extension
20	#WIDTHTHRESHOLD	Width Threshold
21	#PTRESULT	Performance Test Result
22	#PTDATE	Performance Test Date

## Data format group

The *Data format group* contains keywords that describe the instrument. This group is collapsed by default.



Sort	Keyword	Description
	FCS Standard	FCS Version
1	\$MODE	File Mode
2	\$DATATYPE	Data Type
3	\$BYTEORD	Byte Order
4	\$BEGINSTEXT	Begin Text
5	\$ENDSTEXT	End Text
6	\$BEGINDATA	Begin Data
7	\$ENDDATA	End Data

(continued)

Sort	Keyword	Description
8	\$BEGINANALYSIS	Begin Analysis
9	\$ENDANALYSIS	End Analysis
10	\$NEXTDATA	Next Data
6	\$BEGINDATA	Begin Data

## Other group

The *Other group* contains all other keywords not captured within the other groups. This group is collapsed by default.

## Filter configuration (FC) module

### Overview

The *Filter Configuration (FC) module* of the Attune™ Cytometric Software consists of the *File* (“File tab” on page 72) and *Configuration ribbon tab* (“Configuration ribbon tab” on page 513), and the *Filter Configuration View* (“Filter configuration view” on page 526).

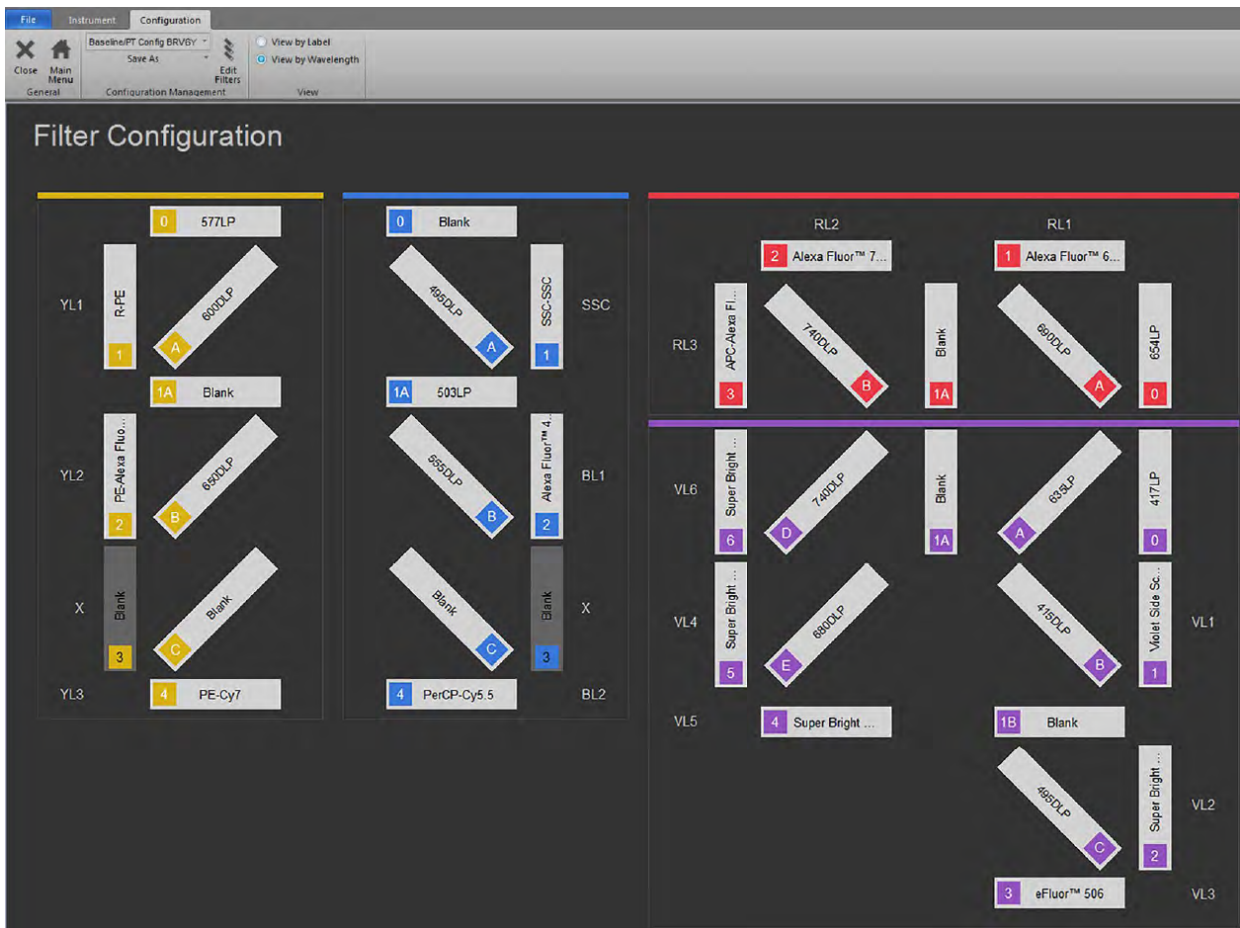


Figure 106 Attune™ NxT BRV6Y configuration

The FC module is used to:

- Select view options for the filters (“View group” on page 524)
- Manage filter labels (“Manage filter labels” on page 524)
- Create, edit, and delete filters (“Edit filters” on page 518)
- Select instrument configuration files (“Select configuration” on page 514)
- Save, load, export, and delete instrument configuration files (“Save as” on page 515)

## Launch the FC module

On the *Instrument ribbon tab* (“Instrument tab” on page 88), click the **Configuration** button to launch the *Filter Configuration (FC) module*.

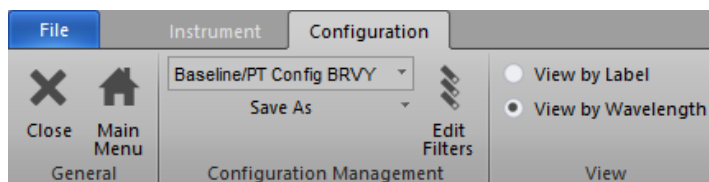


The Configuration button is always visible and enabled except during acquisition.

- When the FC module is opened, all other panels are hidden. The FC module is not resizable or floatable.
- The FC module contains only the File (“File tab” on page 72), Configuration (“Configuration ribbon tab” on page 513), and Instrument (“Instrument tab” on page 88) ribbon tabs. All other ribbons tabs are hidden.

## Configuration ribbon tab

The *Configuration ribbon tab* allows you to manage the filter configuration and to select view options for the filters. It is visible only when the FC module is open.



The Configuration ribbon tab is organized into three functional groups:

- *General* (“Configuration ribbon tab” on page 513)
- *Configuration Management* (“Configuration management group” on page 514)
- *View* (“View group” on page 524)

## General group

*General* group allows you to close the IC module and to return to the Main Menu.

**✕ Close:** Closes the IC module and returns to the previous view before the Instrument Configuration panel was opened. If you have unsaved configuration changes, the *Save As* dialog (“Save as” on page 515) appears.

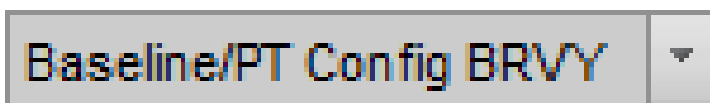
**🏠 Main Menu:** Closes the IC module and returns to the Main Menu. If you have unsaved configuration changes, the *Save As* dialog (“Save as” on page 515) appears.

## Configuration management group

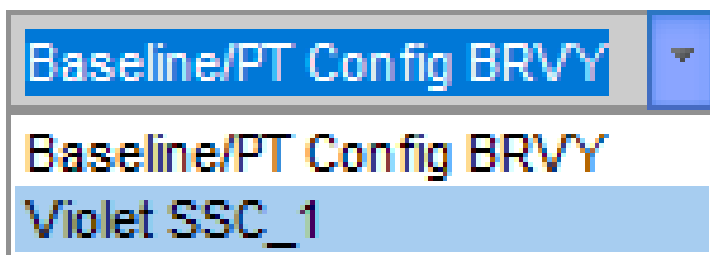
*Configuration Management* group allows you to select instrument configuration, to save, load, export, and delete filter configuration files, and edit the table of filters to be saved as a user-defined configuration.

### Select configuration

**Select configuration** dropdown menu displays the currently selected instrument configuration.



- The default filter configuration is the baseline configuration provided by the instrument manufacturer and is called **Baseline/PT Config <CCC#C>**, where <C> is the first letter of each laser color (**B**lue, **R**ed, **V**iolet, or **Y**ellow) and # corresponds to 4 or 6 channels detected off the Violet laser. An **X** for laser color indicates that one of the lasers is not present.
- The Baseline/PT Configuration is set by the service engineer at time of instrument installation.
- The file containing the default filter configuration cannot be overwritten or deleted.
- To change the filter configuration, select the file for the configuration of interest from the **Select configuration** dropdown menu. The Filter Configuration View (“Filter configuration view” on page 526) is automatically updated to display the filter information for the selected configuration.



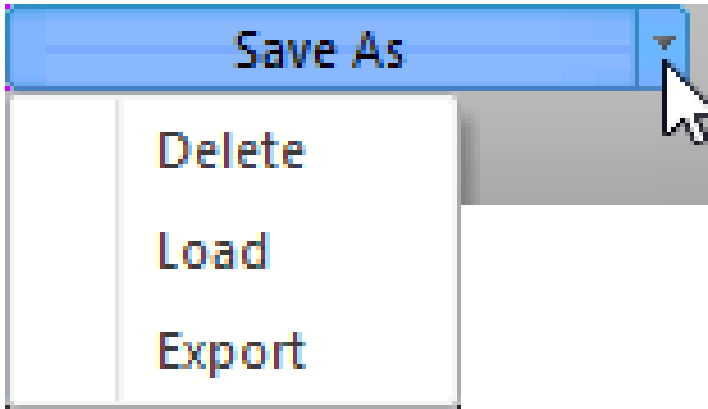
Only filter configurations that match the **Hardware/Virtual Laser Configuration** as defined in the *Options dialog* (“Hardware/Virtual laser configuration” on page 538) are displayed.

- New configurations are user-account specific and can be shared with other users by saving the configuration file, then exporting it to a shared location for import into another user’s account.



## Save as

The Save As button is a split button used for performing the **Save As**, **Load**, **Export**, and **Delete** functions.



**Save As:** Allows you to save the changes to a new file, or overwrite an existing file, if you made changes to the current filter configuration file.

To save changes to a filter configuration file:

1. Click the main portion of the **Save As** button to open the Save As dialog.



2. To save the changes to:
  - A new file – Enter a file name in the text field.
  - An existing file – Select the file from the dropdown menu or enter the name of the existing file in the text field. You can enter up to 50 characters; the following characters are not permitted: \ / : \* < > | ?.

---

**Note:** If you select the default filter configuration file, the **Save** button is disabled.

---

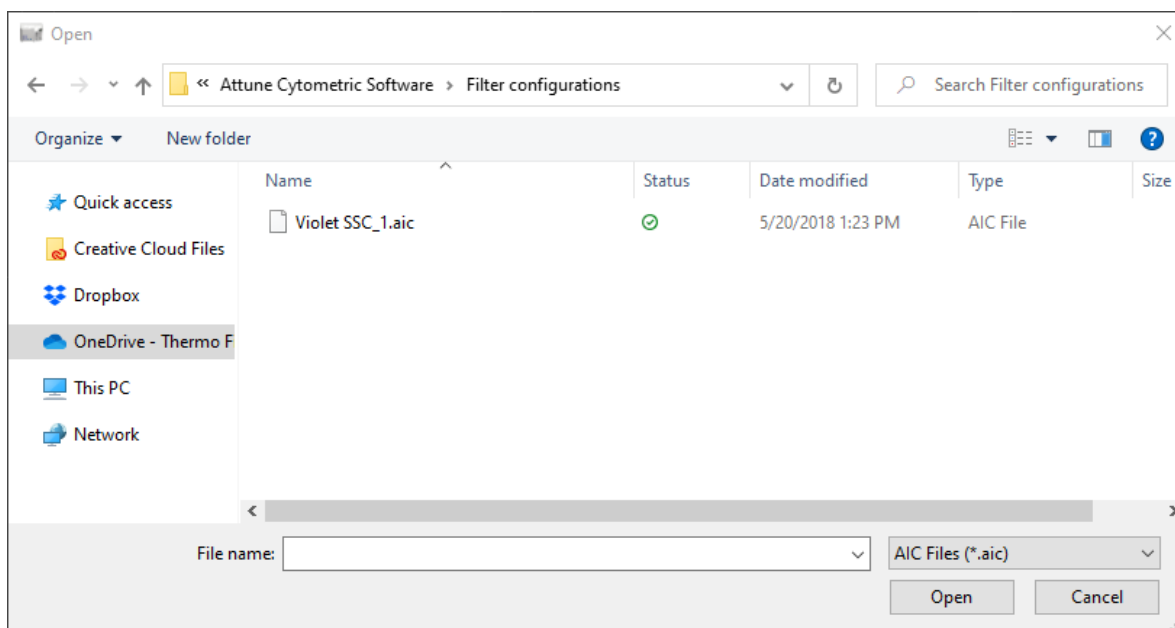
3. Click **Save** (for a new file) or click **Overwrite** (for an existing file) to save the changes.



Alternatively, click **Discard** or **X** to close the *Save As* dialog without saving the changes.

**Load:** Allows you to import a saved filter configuration. To load (import) a filter configuration file:

1. Click the **arrow** next to the **Save As** button, then select **Load**. The *Open* dialog opens. The default location is the location of the last viewed folder.

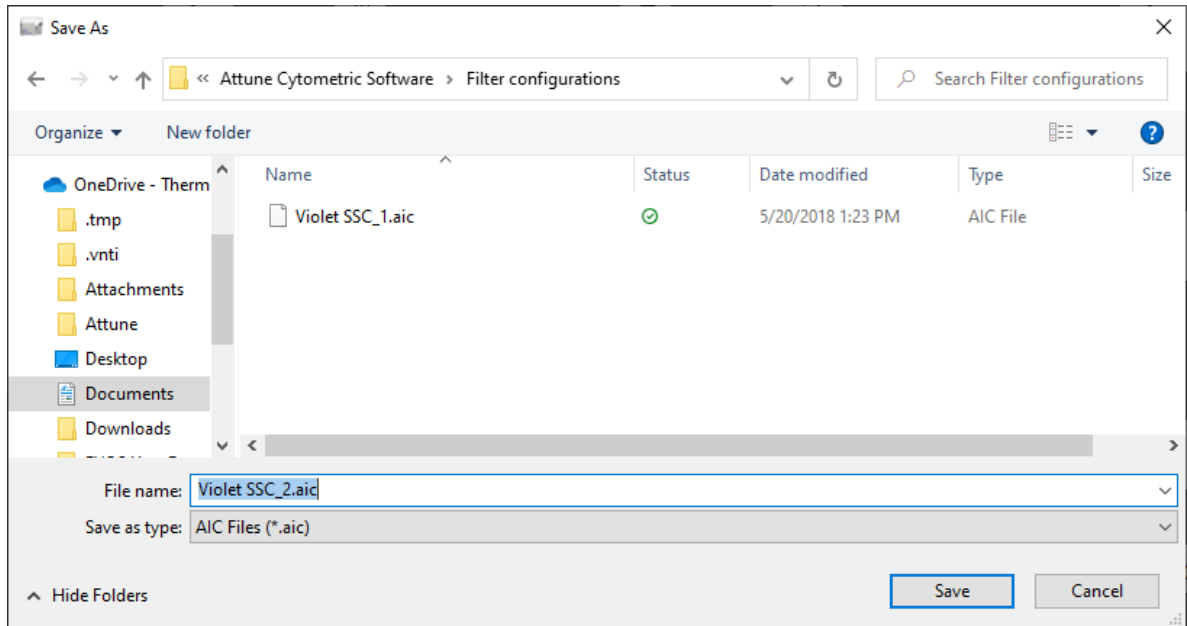


2. Select a file and then click **Import**.

If you select a file that has the same name as a file in your own user list, the software appends an integer onto the file name upon loading.

**Export:** Allows you to export a saved filter configuration file. To export a filter configuration file:

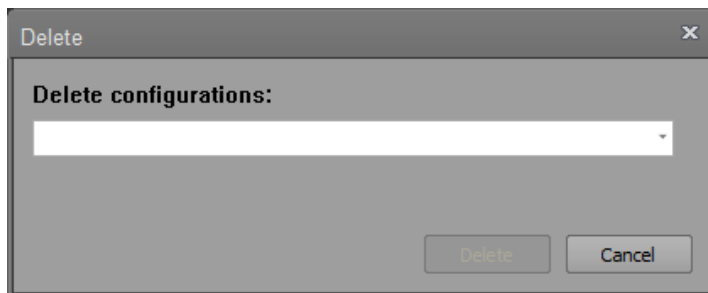
1. Click the **arrow** next to the **Save As** button, then select **Export** to open the *Save As* dialog. The *Save As* dialog automatically opens the location where you last saved files.



2. Browse to a save location, enter a file name, select the **\*.ais** file type, then click **Save**.

**Delete:** Allows you to delete a filter configuration file. To delete a filter configuration file:

1. Click the **arrow** next to the **Save As** button, then select **Delete** to open the *Delete* dialog.



2. By default, the currently loaded filter configuration file is selected. Accept the default or select another file from the dropdown menu.

The dropdown menu displays all available filter configuration files, except for the default filter configuration file.

3. Click **Delete** to delete the selected file and close the dialog box.

Alternatively, click **Cancel** or **X** to close the *Delete* dialog without deleting the file.

---

**Note:** If you delete the currently loaded filter configuration file, the software updates the view to the default filter configuration file.

---

## Edit filters

**Edit Filters** button opens the *Edit Filters dialog* (“Edit filters dialog” on page 519). The dialog contains the *Filters table*, the contents of which are based on the filter configuration selected from the **Select configuration** dropdown menu (“Configuration ribbon tab” on page 513).



Using the Edit Filters dialog, you can:

- Define a new filter (“Define a new filter” on page 521)
- Edit an existing filter (“Edit an existing filter” on page 523)
- Delete an existing filter (“Delete a filter” on page 523)
- Manage filter labels (“Manage filter labels” on page 524)

---

**Note:** Only *Advanced™ Users* and *Administrators* (“Account permissions” on page 44) can perform the tasks in the *Edit Filters* dialog.

---

## Edit filters dialog

Using the tools available in the Edit Filters dialog, authorized users can define new filter sets, change the filters that can be applied to each instrument channel, and manage filter labels.

**Filters**

Laser Line	Filter Type	Wavelength	Primary Label	Delete
Blue	Collection	488/10	SSC-SSC	N/A
Blue	Directing	495DLP		N/A
Blue	Directing	503LP		N/A
Blue	Collection	530/30	FITC	N/A
Blue	Directing	555DLP		N/A
Blue	Collection	590/40	PE-Texas Red™	N/A
Blue	Directing	650DLP		N/A
Blue	Collection	695/40	PerCP	N/A
Blue	Directing	Blank		N/A
Blue	Directing	Blank		N/A
Blue	Directing	Blank		N/A
Red	Directing	654LP		N/A
Red	Collection	670/14	APC	N/A
Red	Directing	690DLP		N/A
Red	Collection	720/30	Alexa Fluor™ 700	N/A
Red	Directing	740DLP		N/A
Red	Collection	780/60	APC-Alexa Fluor™ 750	N/A
Red	Directing	Blank		N/A
Red	Directing	Blank		N/A

## Filters table

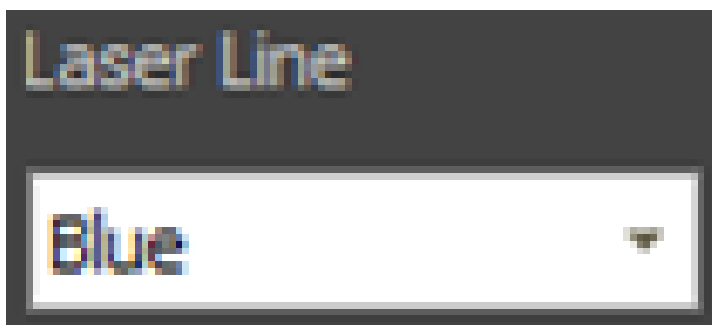
The Filters table displays the filters for the currently selected filter configuration, with each filter taking one row of the table.

- The columns in the table are listed from left to right as follows: Laser Line, Filter Type, Wavelength, Primary Label, and Delete button.
- By default, the filters are sorted by the laser line group, and the secondary sort order is wavelength.
- Authorized users can edit the contents of this table as described below. Changes to the table are saved to the user-defined filter configuration.

## Edit filters tools

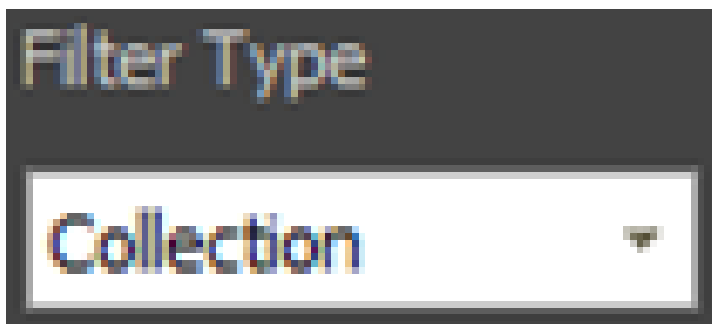
The upper portion of the Edit Filters dialog contains the tools for defining a new filter or editing an existing filter. It contains the following options:

- **Laser Line dropdown:** Allows you to select the laser to which the filter will be assigned. The default is the first laser in the configuration file.

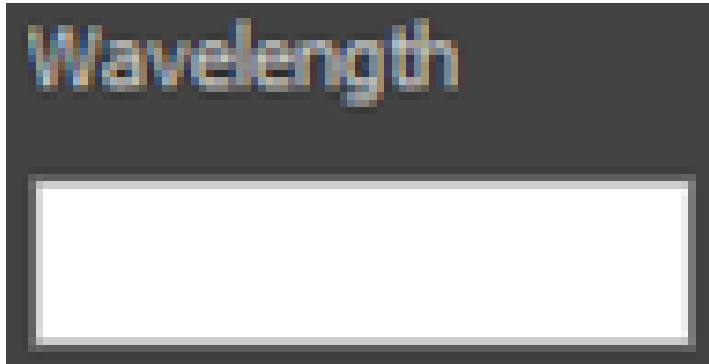


When an Attune™ system is detected or selected within the virtual mode, only the lasers which are installed in that instrument are available.

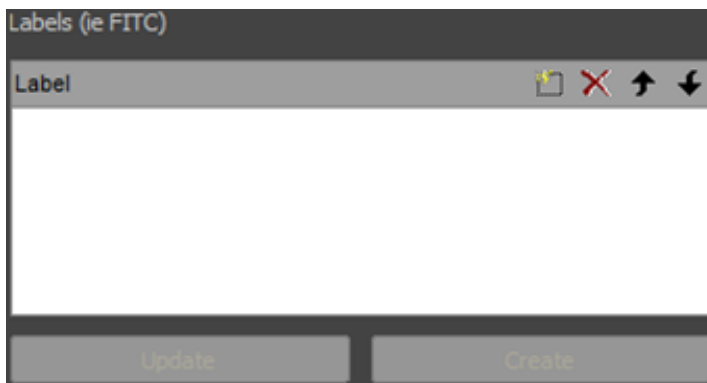
- **Filter Type dropdown:** Allows you to select the filter type. You can choose between **Band Pass**, **Long Pass**, **Short Pass**, and **Dichroic**. By default, Band Pass is selected.



- **Wavelength:** Allows you to type in the filter specifications. You can enter up to 10 characters to the text field. By default, this field is blank.



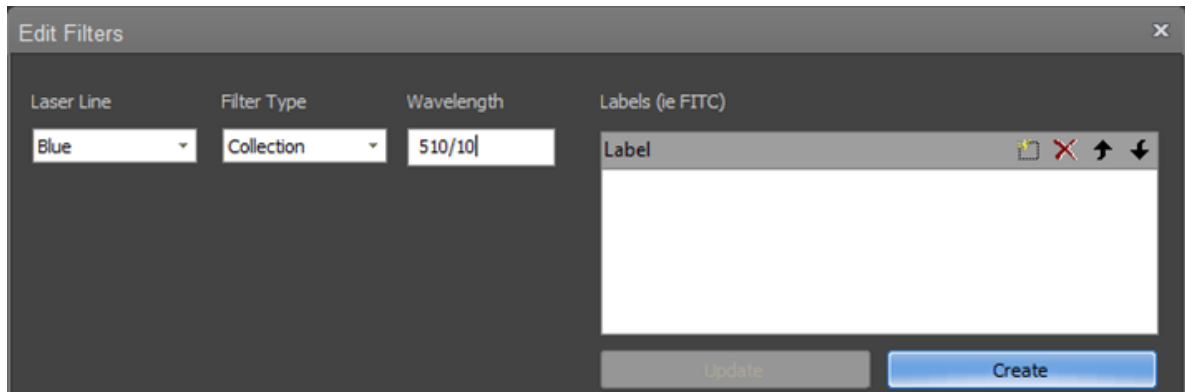
- **Labels table:** Allows you to assign and manage filter labels using the tools available in the table header (see “Manage filter labels” on page 524). You can enter up to 50 characters per line to save as a label for the selected filter. By default, this table is blank.  
**Update** and **Create** buttons are contextual; they are enabled only when the appropriate action has been taken in the Edit Filters dialog.



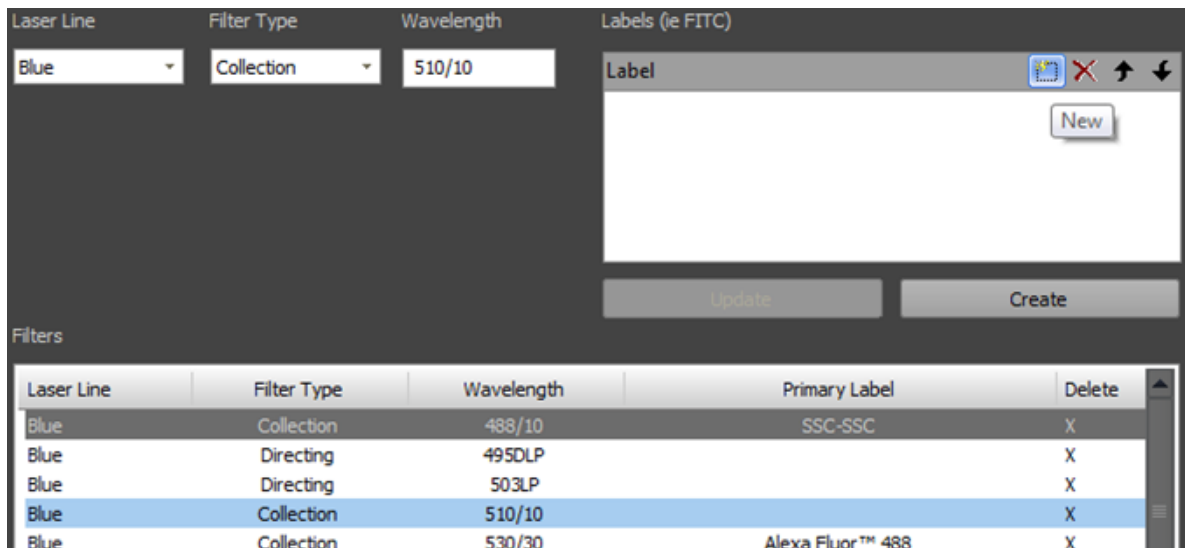
### Define a new filter

1. Click **Edit Filters** to open the *Edit Filters* dialog.
2. In the **Laser Line** dropdown menu, select the laser to which you want to assign the new filter.
3. In the **Filter Type** dropdown menu, select the desired filter type.

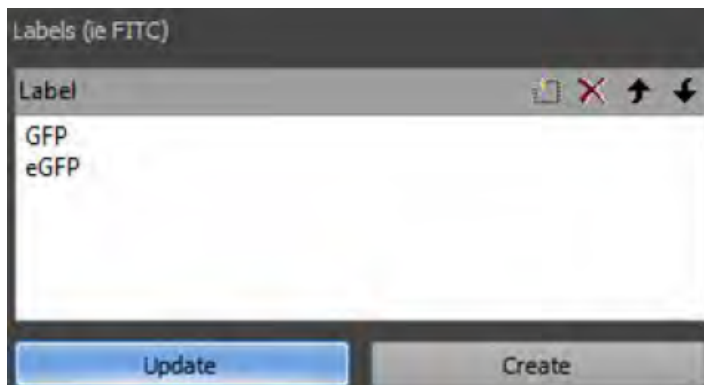
- In the **Wavelength** field, enter the filter specifications. You can enter up to 10 characters.



- Click **Create**. The software adds the new filter to the **Filters** table.
- To assign a *primary label* (i.e., the default display name) for the filter, select the new filter from the Filters table, then click **New** in the Labels table.



- Type the name for the new filter in the Labels table, then click **Update**. You can create additional labels by repeating this process.





8. If needed, you can edit the label, or create a new label (see “Manage filter labels” on page 524).
9. Click **OK** to apply the changes and close the *Edit Filters* dialog. The software saves the changes to the currently selected instrument configuration file. Changes will be lost if the **OK** button is not clicked before closing the menu.  
Alternatively, click **Cancel** to close the *Edit Filters* dialog without applying the changes.

### Edit an existing filter

1. Click **Edit Filters** to open the *Edit Filters* dialog.
2. In the **Filters** table, select the filter to edit. The software populates the fields with the selected filter’s information.
3. For user-defined filters, you can make the following changes:
  - In the **Laser Line** dropdown menu, select the laser to assign the filter to.
  - In the **Filter Type** dropdown menu, select the desired filter type.
  - In the **Wavelength** field, enter the filter specifications. You can enter up to 10 characters.

---

**Note:** The Laser Line, Filter Type, and Wavelength fields are disabled for the filters in the system default configuration.

---

4. Assign the primary label for the selected filter by selecting the label of interest in the **Labels** table, and then using the up arrow in the header to move it to the top of the list.  
If needed, you can edit the label, or create a new label (see “Manage filter labels” on page 524).
5. Click **Update**.
6. Click **OK** to apply the changes and close the *Edit Filters* dialog. The software saves the changes to the currently selected instrument configuration file. Changes will be lost if the **OK** button is not clicked before closing the menu.  
To close the *Edit Filters* dialog without applying the changes, click **Cancel**.

### Delete a filter

You can only delete filters that were created by a user; filters installed with the system cannot be deleted.

1. Click **Edit Filters** to open the *Edit Filters* dialog.
2. In the **Filters** table, select the filter to delete and click **X** in the Delete column.
3. Click **OK** to apply the changes and close the *Edit Filters* dialog. The software saves the changes to the currently selected instrument configuration file. Changes will be lost if the **OK** button is not clicked before closing the menu.  
To close the *Edit Filters* dialog without applying the changes, click **Cancel**.

## Manage filter labels

You can arrange, create, edit, and delete filter labels in the **Labels** table.

1. Click **Edit Filters** to open the *Edit Filters* dialog.
2. In the **Filters** table, select the filter to edit. The software populates the fields with the selected filter's information.
3. In the **Labels** table, perform the following tasks as needed:
  - Click on a label to select it. To select multiple labels at a time, press **Ctrl** or **Ctrl+Shift** when selecting the labels.
  - To arrange labels in the table, select a label, then use the up or down arrows in the header (⬆️ ⬇️) to move the selected label.  
The label at the top of the list is considered the *primary label* and will be used as the default display name for the selected filter.
  - To create a new label, click **Add** (➕) in the header or double-click the white space within the table, and then enter a label name. You can enter up to 50 characters; the following characters are not permitted: \ / : \* < > | ?.
  - To edit an existing label, double-click the label and enter changes.
  - To delete a label, select it, and then click **Delete** (✖️) in the header.

## View group

View group is used for displaying the filters in the filter configuration by primary label or by assigned wavelength. It consists of **View by label** and **View by wavelength** radio buttons. Selecting one option deselects the other.



- The default selection is View by wavelength.
- The selection of View by label or View by wavelength is remembered as a user setting and persists on a per user basis.

## Default filter label list

The following tables list the filter labels that are displayed for each channel for the available system default configurations. The naming convention for the system default filter configuration is as follows:

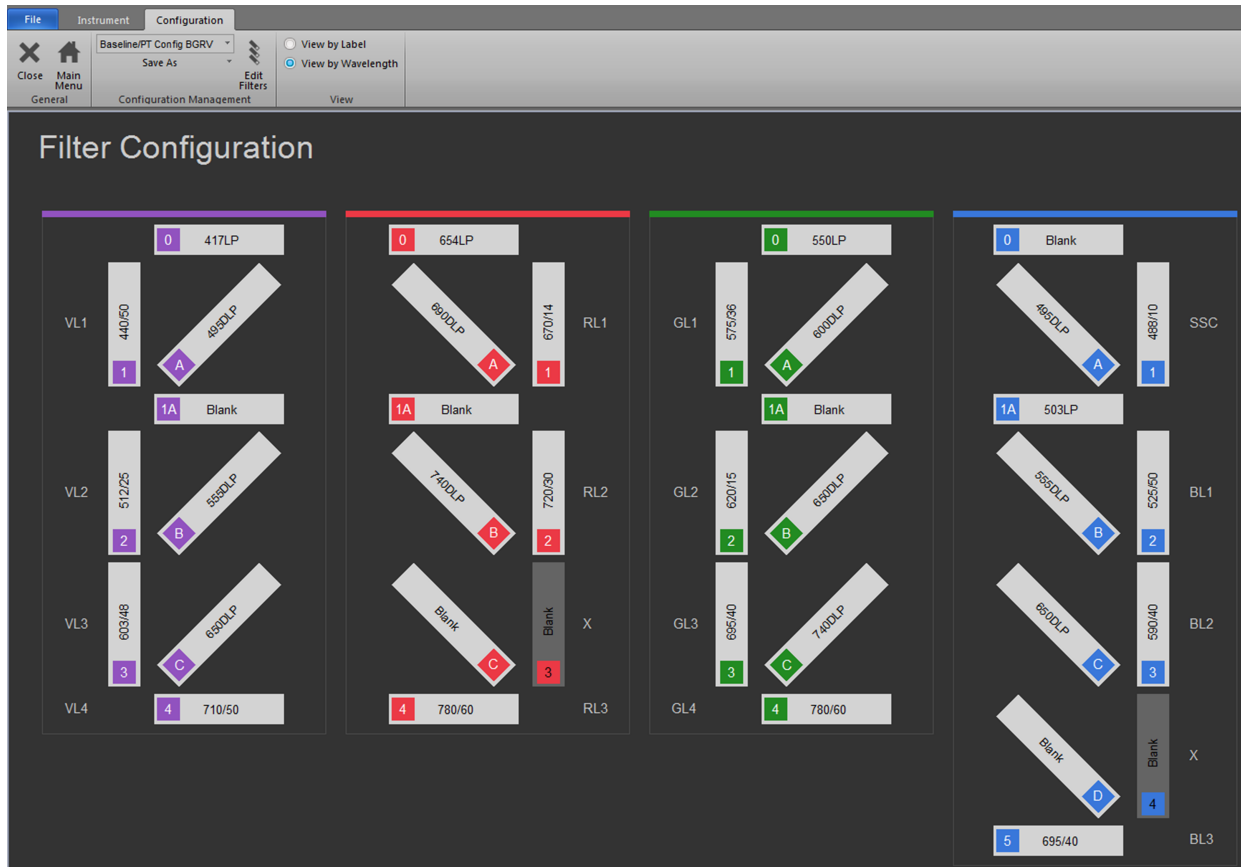
**Baseline/PT Config <CCCC#>**, where **<C>** is the first letter of each laser color (**B**lue, **G**reen, **Y**ellow, **R**ed, or **V**iolet) and **#** corresponds to 4 or 6 channels detected off the Violet laser. An **X** for laser color indicates that one of the lasers is not present.

Configuration	BXXX	BGXX	BYXX	BRXX	BV4XX	BV6XX	BGRX	BRV4X	BGV4X	BYRX	BYV4X	BRV6X	BYRV6	BGRV4	BYRV4	
No. of detectors	4	7	7	7	7	9	10	10	11	11	11	12	14	14	14	
Laser	Channel	Emission filter (nm)														
Blue	BL1	530/30	525/50	530/30	530/30	530/30	530/30	525/50	530/30	525/50	530/30	530/30	530/30	530/30	525/50	530/30
	BL2	574/26	590/40	590/40	574/26	574/26	574/26	590/40	574/26	590/40	574/26	590/40	574/26	695/40	590/40	590/40
	BL3	695/40	695/40	695/40	695/40	695/40	695/40	695/40	695/40	695/40	695/40	695/40	695/40		695/40	695/40
	BL4	780/60			780/60				780/60		780/60					
Green	GL1		575/36					575/36		575/36					575/36	
	GL2		620/15					620/15		620/15					620/15	
	GL3		695/40					695/40		695/40					695/40	
	GL4		780/60					780/60		780/60					780/60	
Yellow	YL1			585/16							585/16	585/16		585/16		585/16
	YL2			620/15							620/15	620/15		620/15		620/15
	YL3			695/40							695/40	695/40		780/60		695/40
	YL4			780/60							780/60	780/60				780/60
Red	RL1				670/14			670/14	670/14		670/14		670/14	670/14	670/14	670/14
	RL2				720/30			720/30	720/30		720/30		720/30	720/30	720/30	720/30
	RL3				780/60			780/60	780/60		780/60		780/60	780/60	780/60	780/60
Violet	VL1					440/50	450/40		440/50	440/50		440/50	450/40	450/40	440/50	440/50
	VL2					512/25	525/50		512/25	512/25		512/25	525/50	525/50	512/25	512/25
	VL3					603/48	610/20		603/48	603/48		603/48	610/20	610/20	603/48	603/48
	VL4					710/50	660/20		710/50	710/50		710/50	660/20	660/20	710/50	710/50
	VL5						710/50						710/50	710/50		
	VL6						780/60						780/60	780/60		

## Filter configuration view

### Overview

The *Filter Configuration View* displays an interactive diagram of the filter array for the current instrument model, which allows you to assign appropriate filters to specific filter locations.



- Only detectors that are part of the active instrument model are displayed.
- The filter array reflects the **Hardware/Virtual Laser Configuration** as defined in the Configuration *Options dialog* (“Hardware/Virtual laser configuration” on page 538) and can be used as a quick reference.
- Clicking one of the filter locations opens a dropdown menu, which contains all unassigned filters for that laser line and type. You can select any of these to assign it to that filter location. Filters marked as **BP** (Band Pass) and **LP** (Long Pass) can be selected at the detector output.
- Filters marked as **Dichroic** are only available in the positions that accept dichroic filters.
- Default Instrument Configuration panels for each laser configuration are shown on “Default instrument configurations” on page 527.

**Note:** Available laser configurations depend on the type of your instrument (Attune™ NxT or the Attune™ CytPix™ Flow Cytometer).

## Default instrument configurations

The following images show the Instrument Configuration panel for each laser configuration that is available for both Attune™ NxT and Attune™ CytPix™ Flow Cytometers (“Available for Attune™ NxT and Attune™ CytPix™ flow cytometers” on page 527), for Attune™ NxT Flow Cytometers only (“Available for the Attune™ NxT flow cytometer only” on page 531), and for Attune™ CytPix™ Flow Cytometers only (“Available for the Attune™ CytPix™ flow cytometer only” on page 535).

### Available for Attune™ NxT and Attune™ CytPix™ flow cytometers

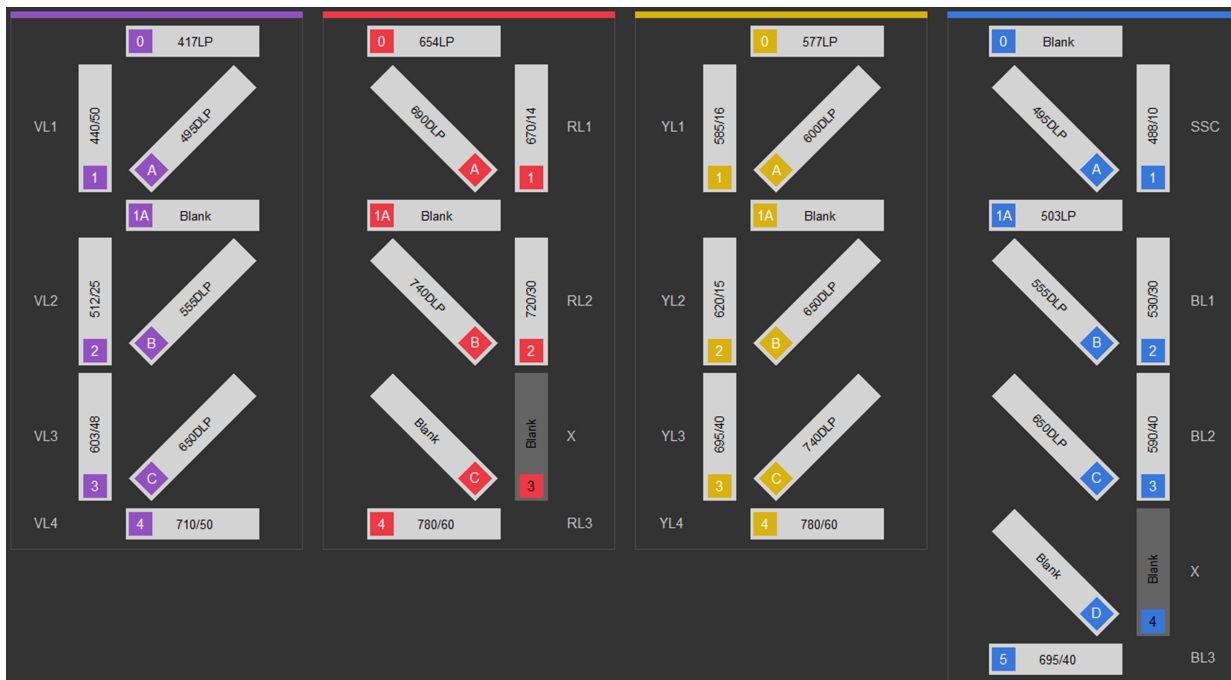


Figure 107 BYRV4 (Blue, Yellow, Red, Violet)

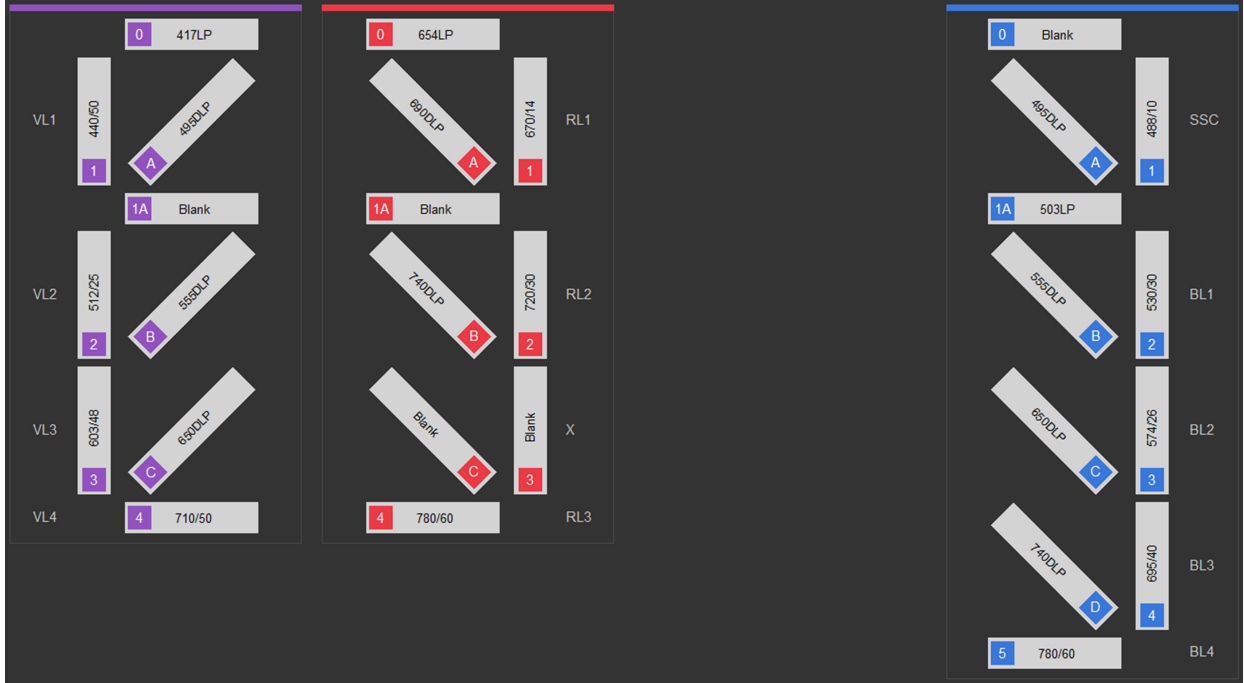


Figure 108 BRV4 (Blue, Red, Violet)



Figure 109 BYV4 (Blue, Yellow, Violet)

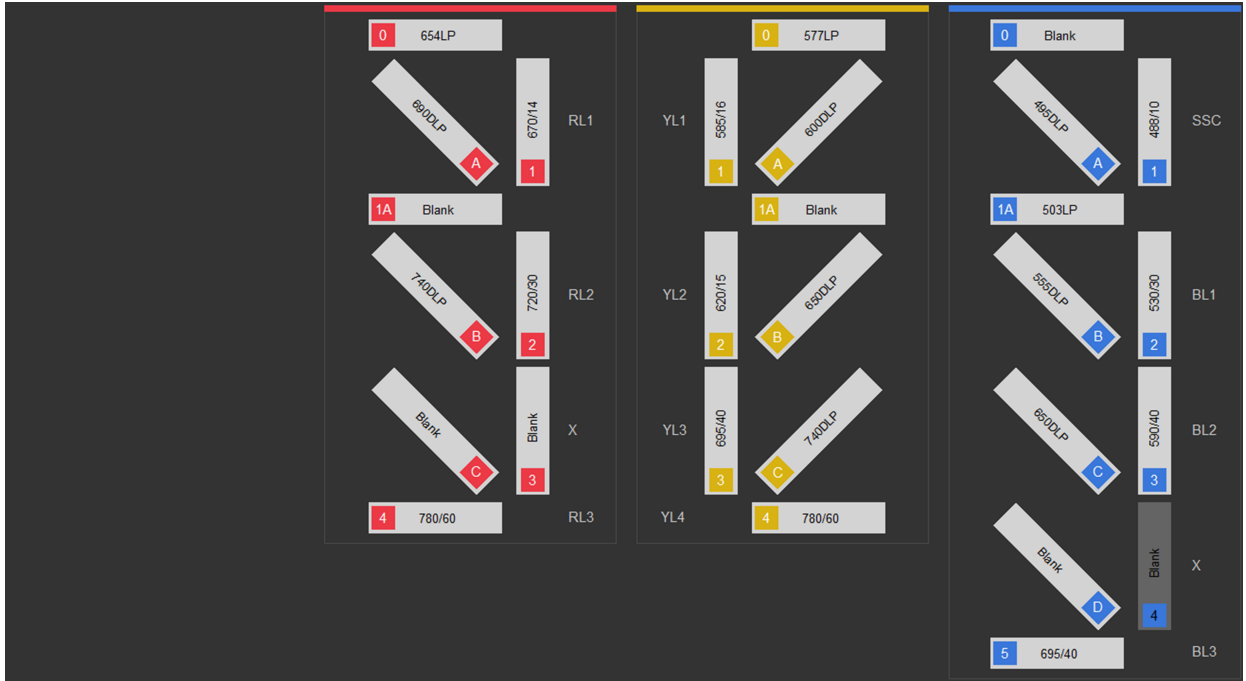


Figure 110 BYR (Blue, Yellow, Red)

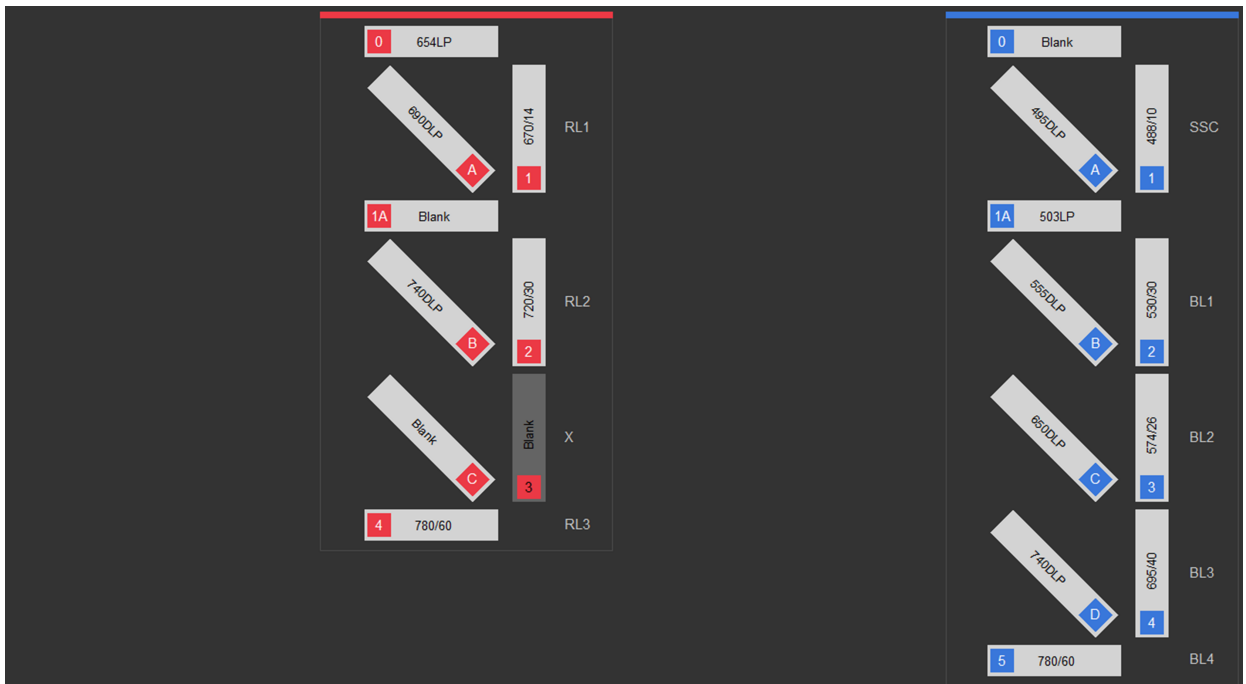


Figure 111 BR (Blue, Red)

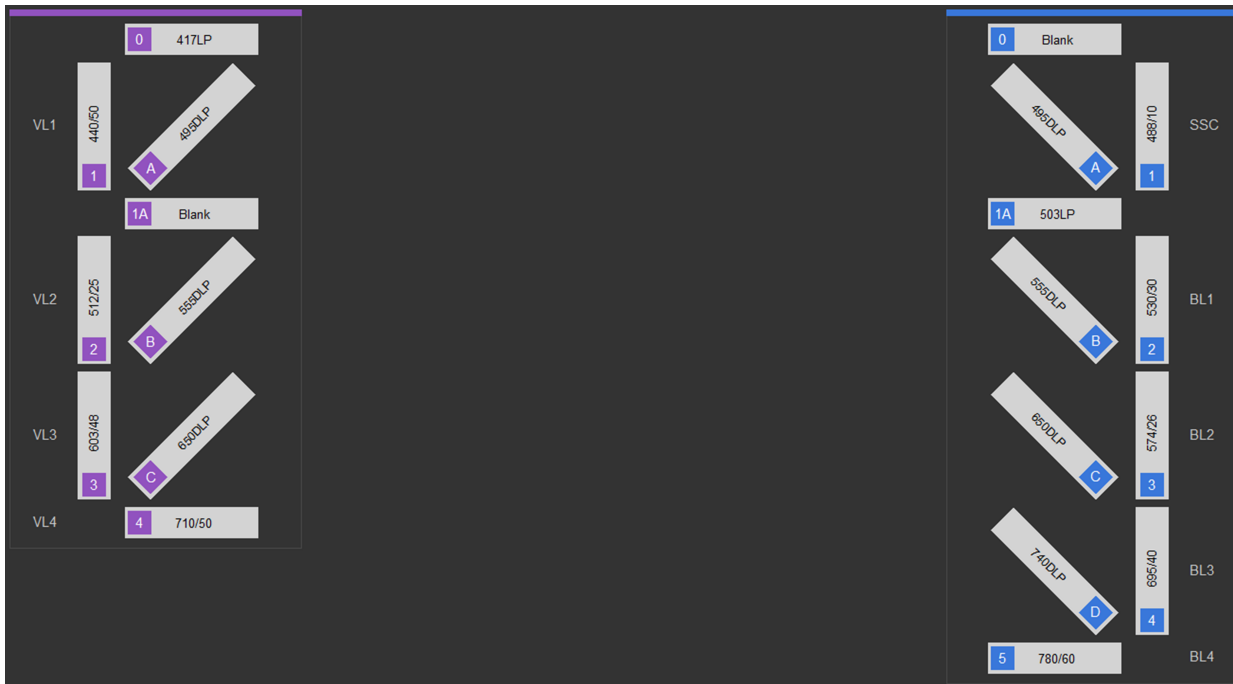


Figure 112 BV4 (Blue, Violet)

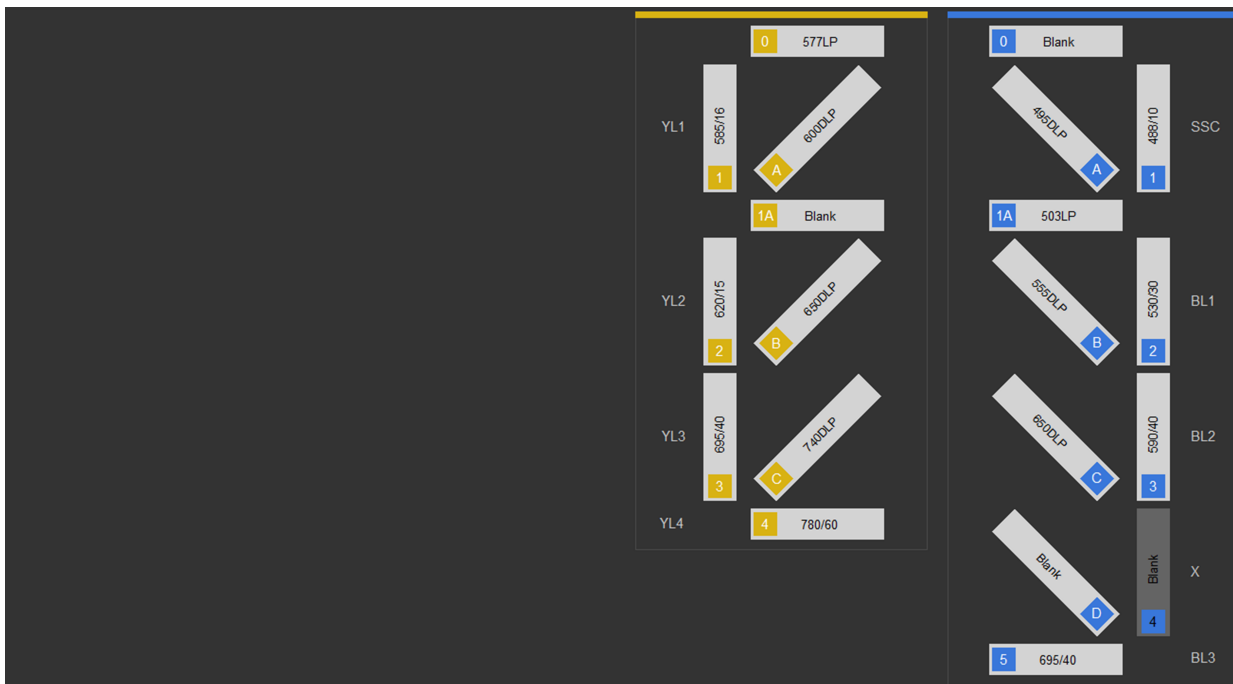


Figure 113 BY (Blue, Yellow)



Available for the Attune™ NxT flow cytometer only

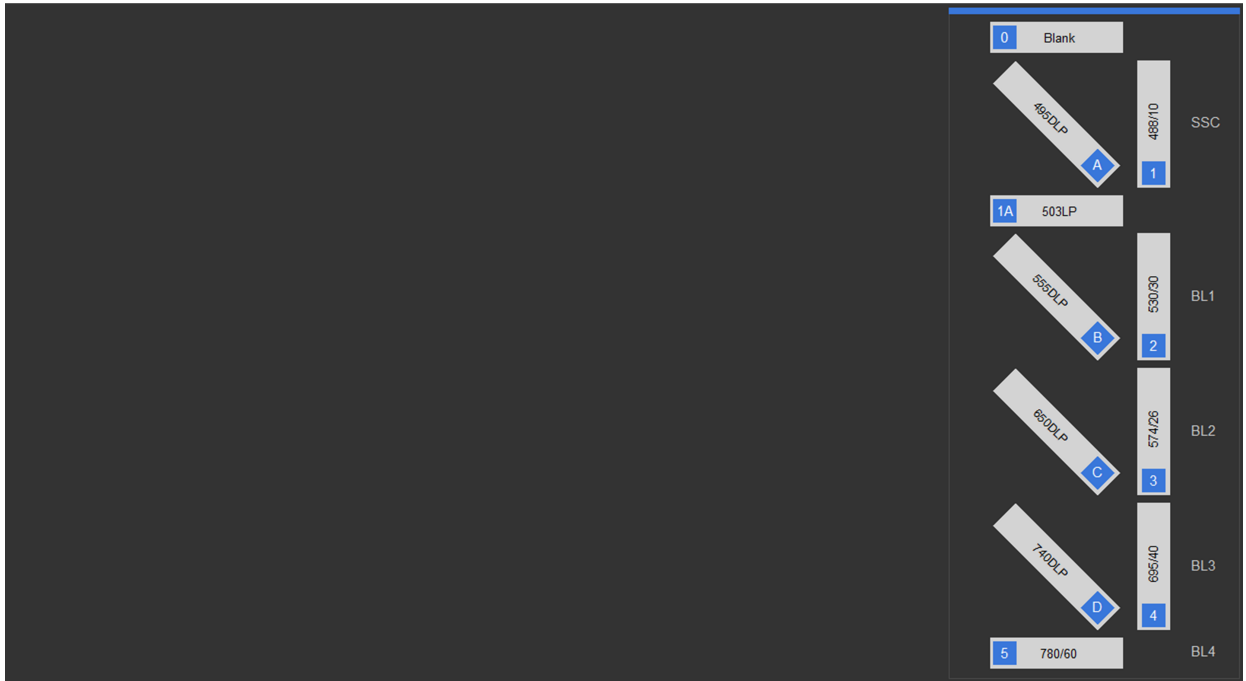


Figure 114 B (Blue)

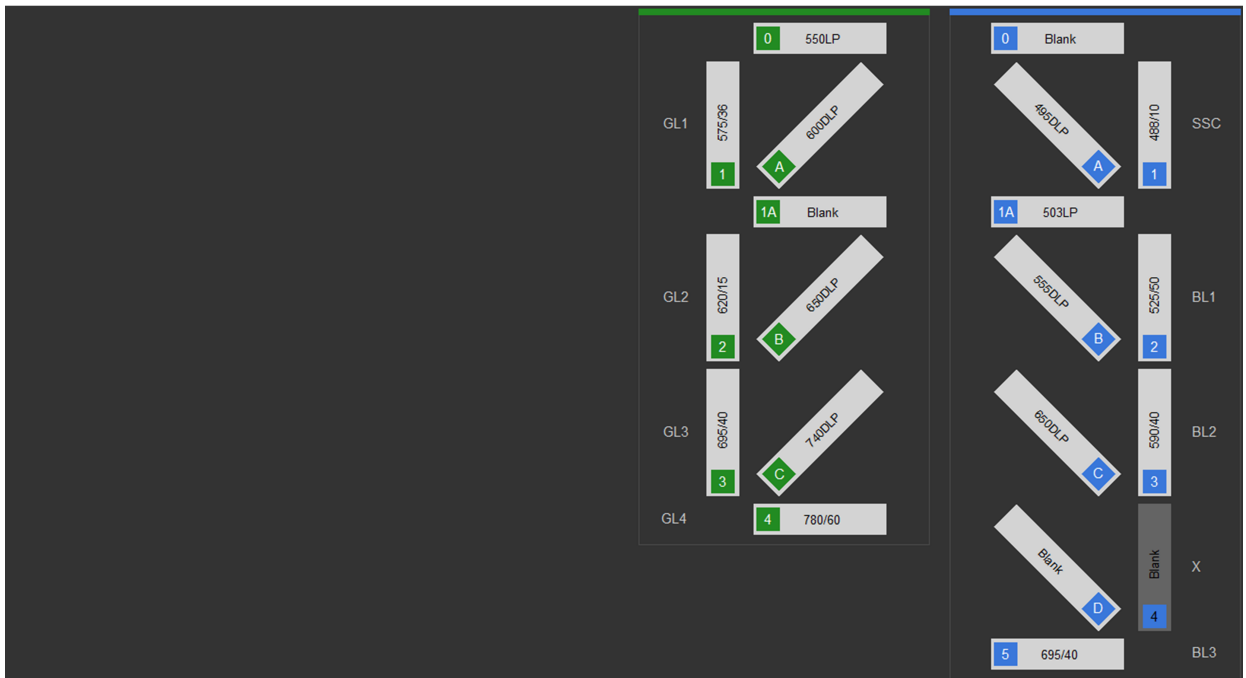


Figure 115 BG (Blue, Green)

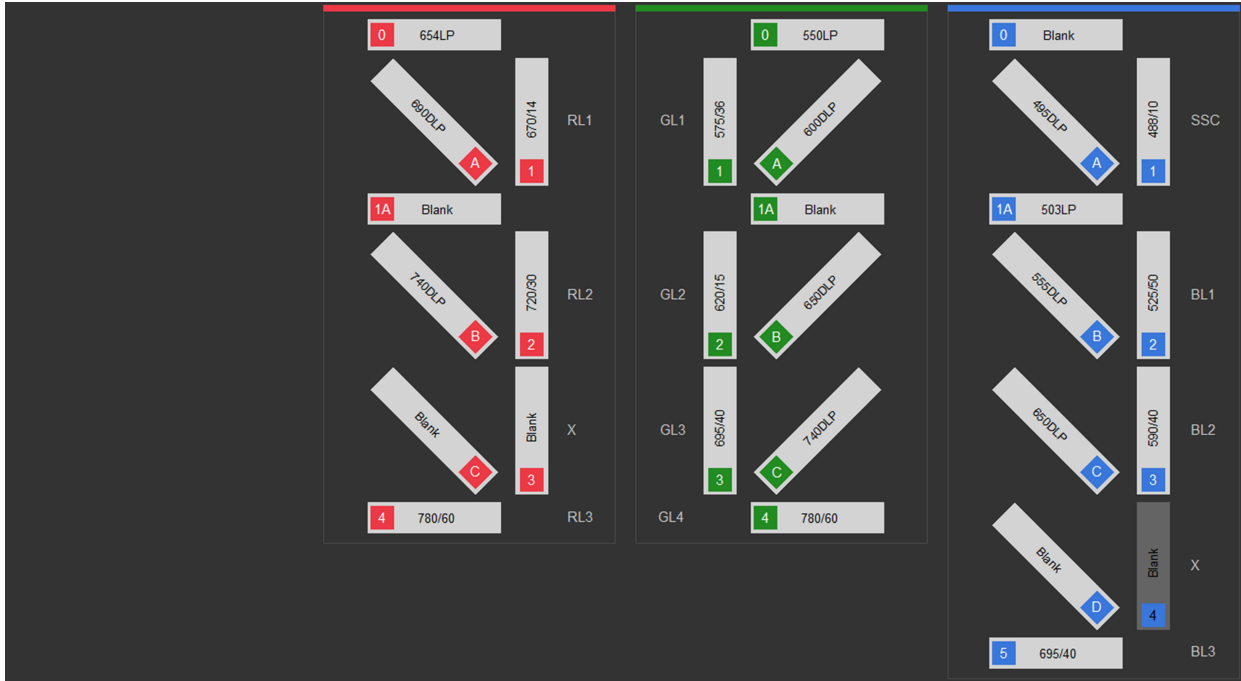


Figure 116 BGR (Blue, Green, Red)

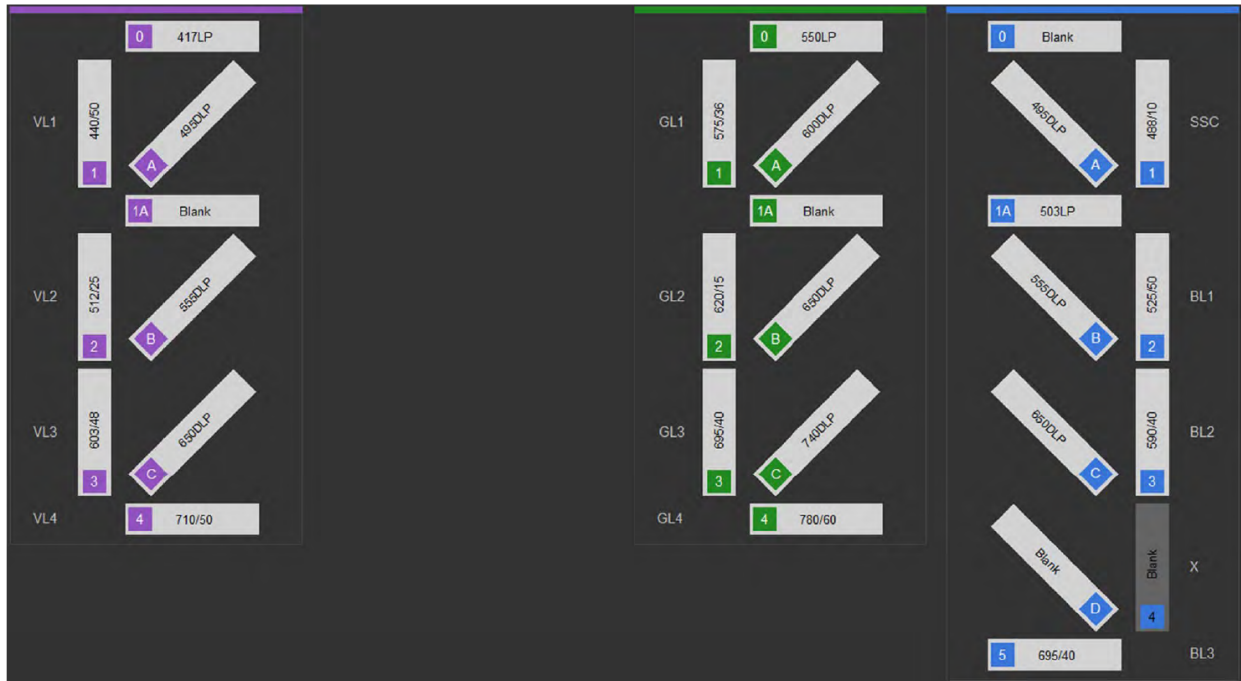


Figure 117 BGV4 (Blue, Green, Violet)

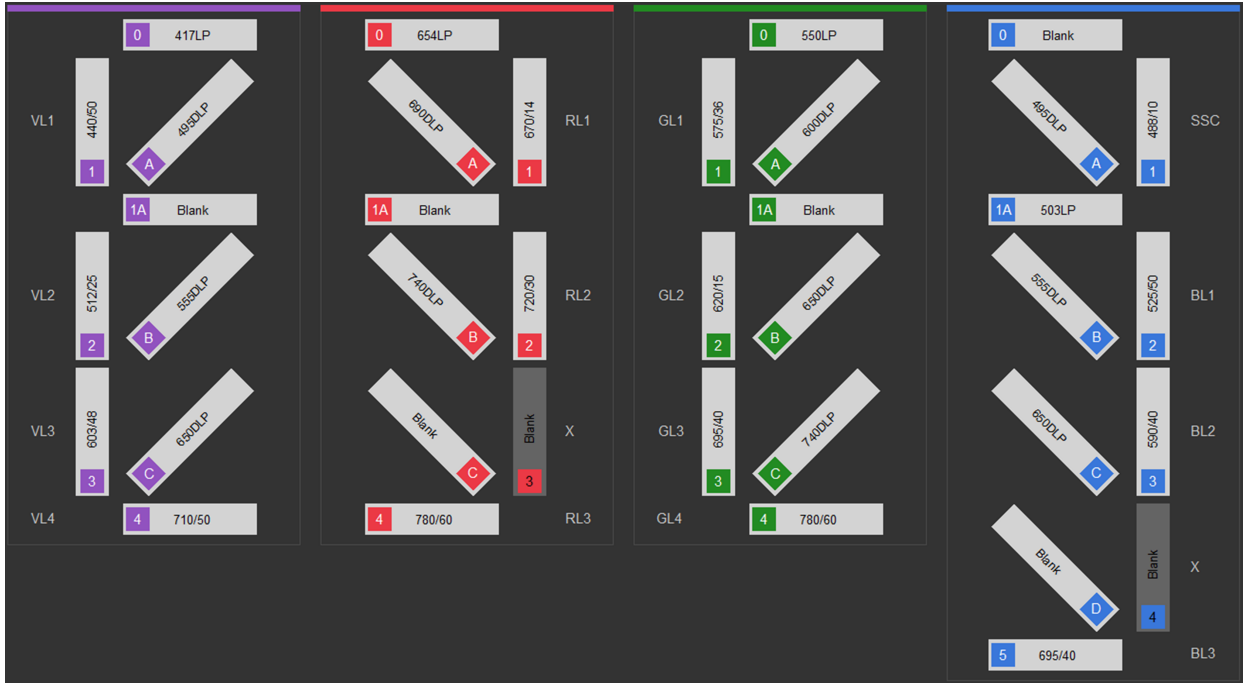


Figure 118 BGRV4 (Blue, Green, Red, Violet)

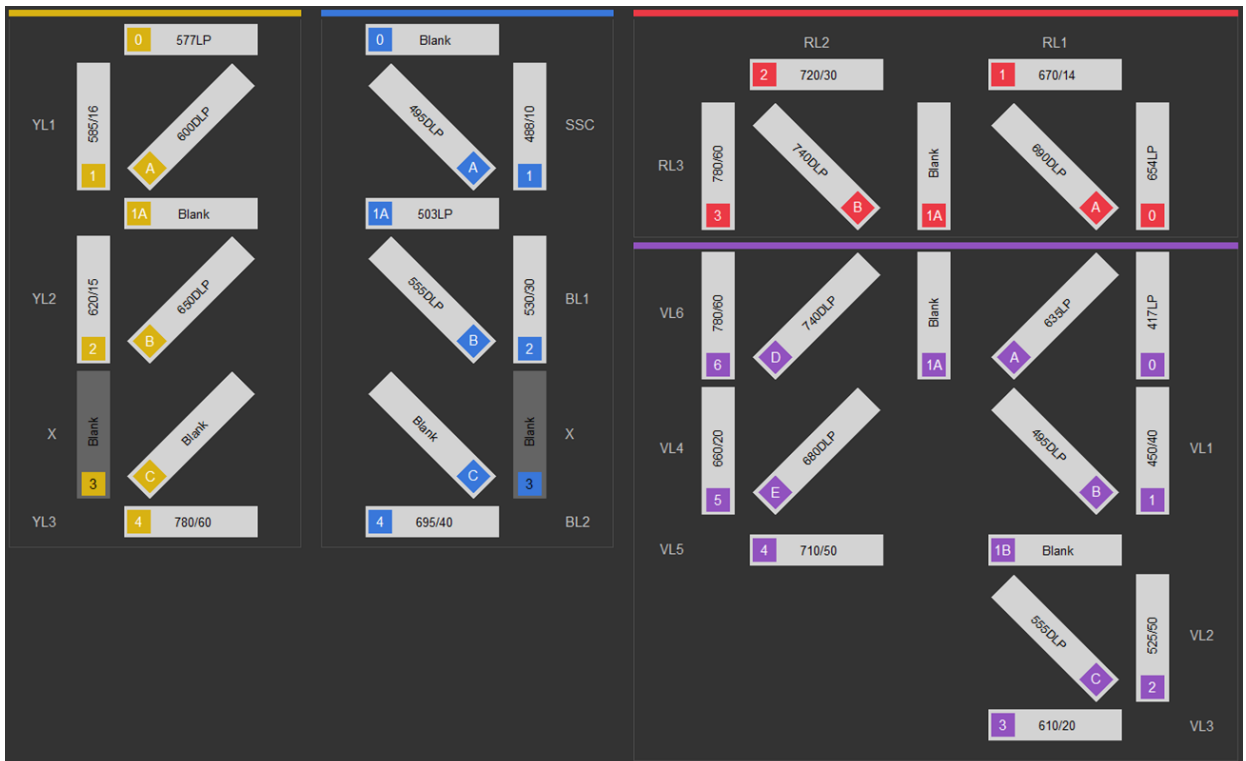


Figure 119 BYRV6 (Blue, Yellow, Red, Violet 6)

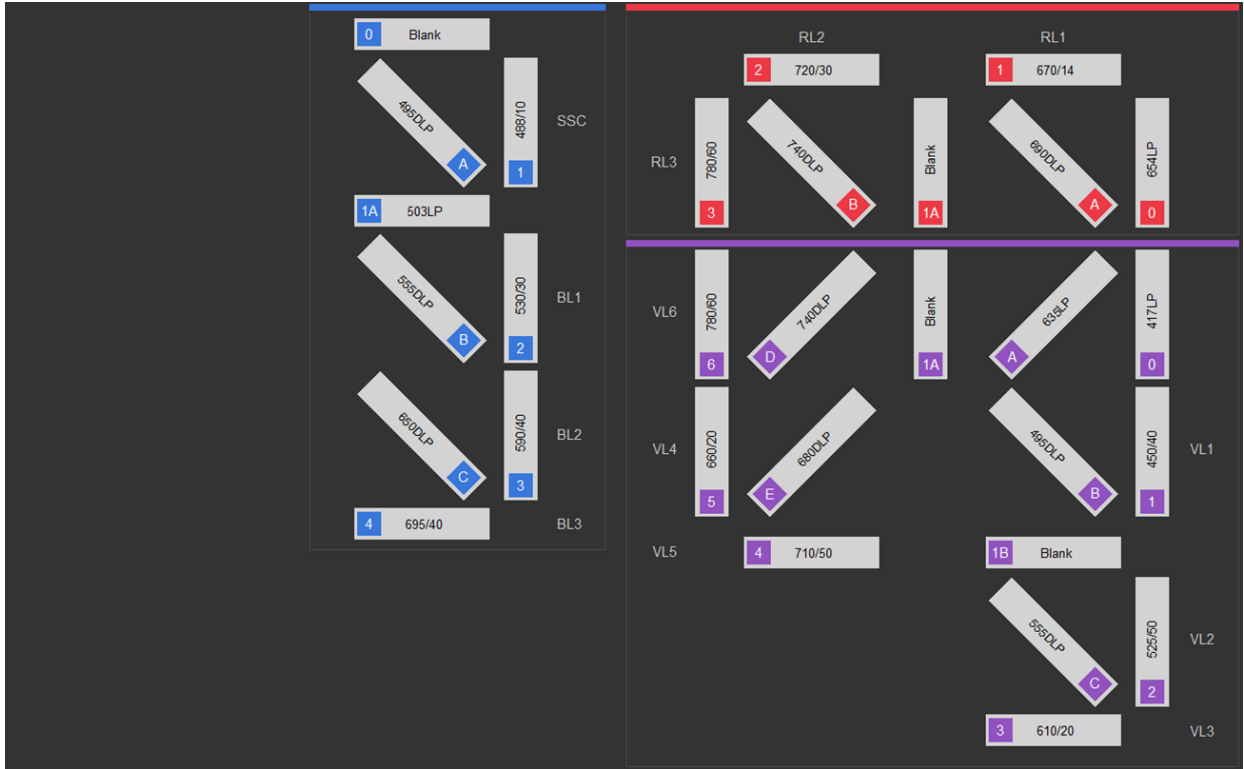


Figure 120 BRV6 (Blue, Red, Violet 6)

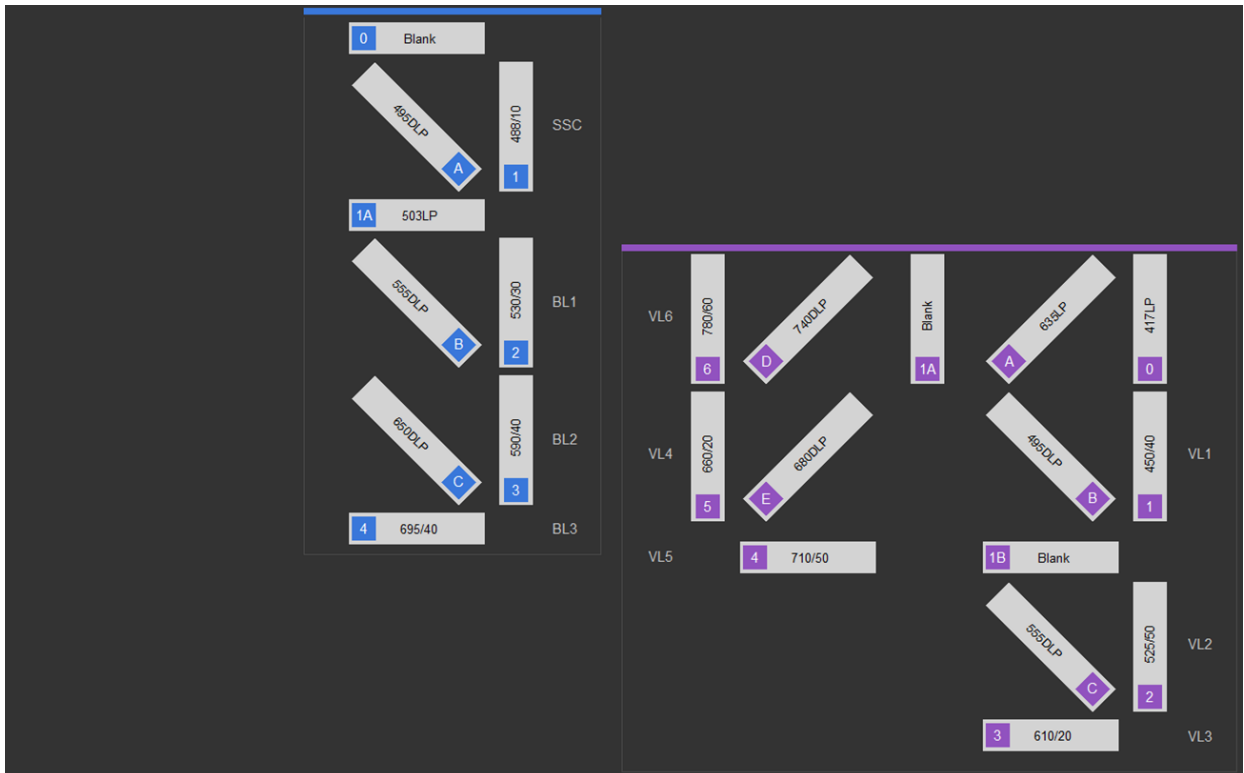


Figure 121 BV6 (Blue, Violet 6)

Available for the Attune™ CytPix™ flow cytometer only

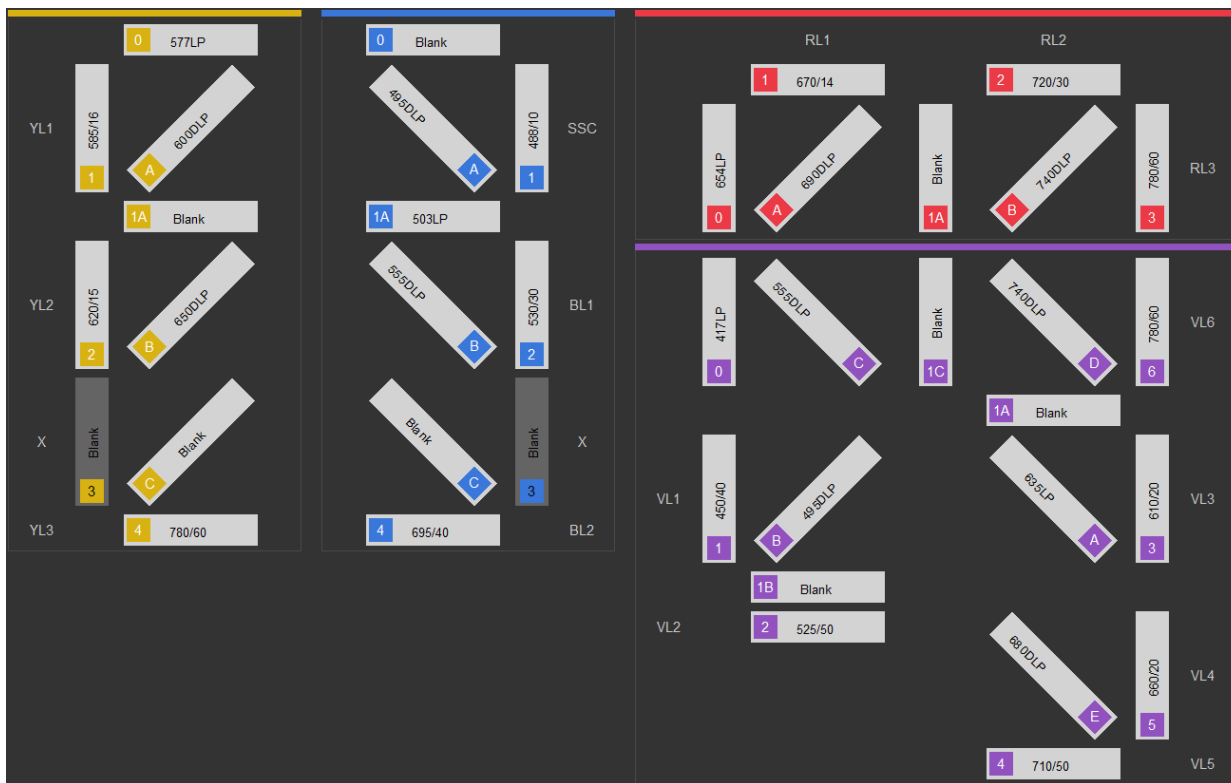


Figure 122 BYRV6 (Blue, Yellow, Red, Violet 6)

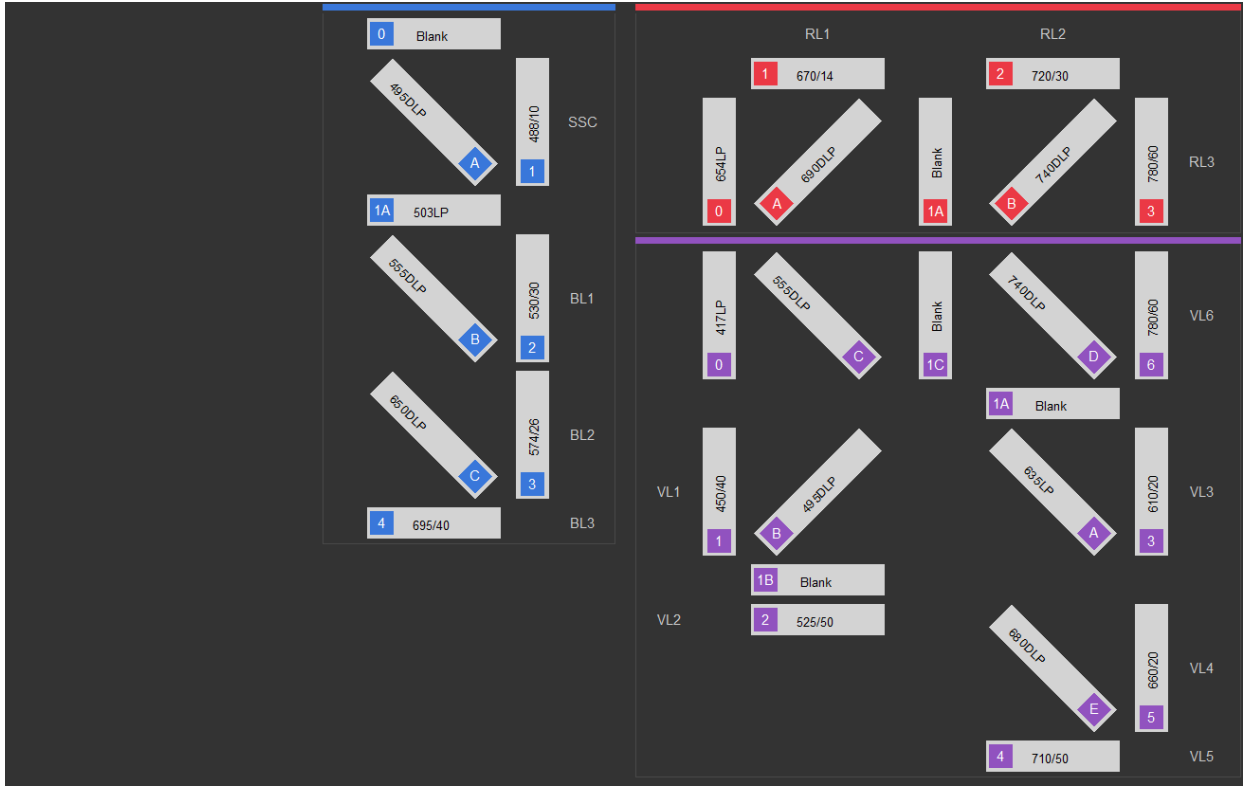


Figure 123 BRV6 (Blue, Red, Violet 6)

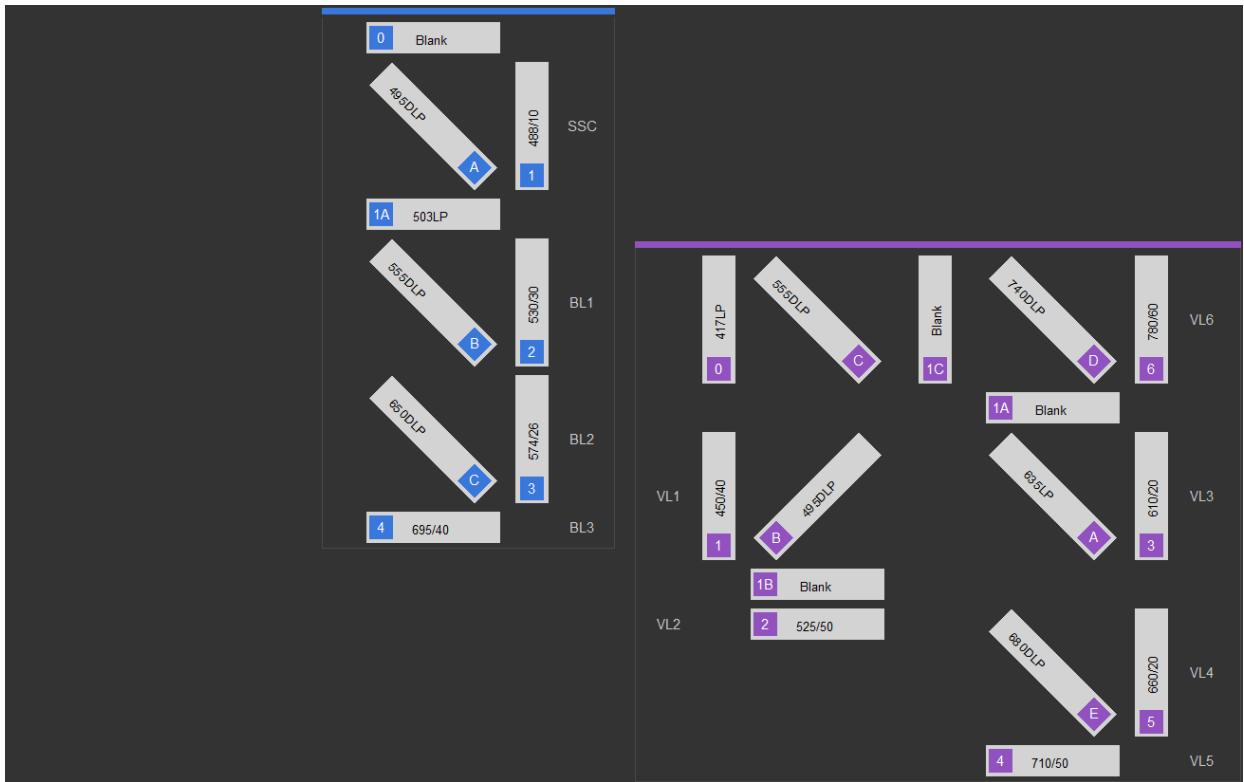


Figure 124 BV6 (Blue, Violet 6)

**Note:** In Attune™ CytPix™ Flow Cytometers, the orientation of the Violet 6 laser configuration is reversed to accommodate the image collection optics.

## Configuration options

### Overview

*Configuration* options allow you to create virtual configuration settings for working in analysis-only conditions, and to manage the bubble sensor and set the sample conservation mode property.

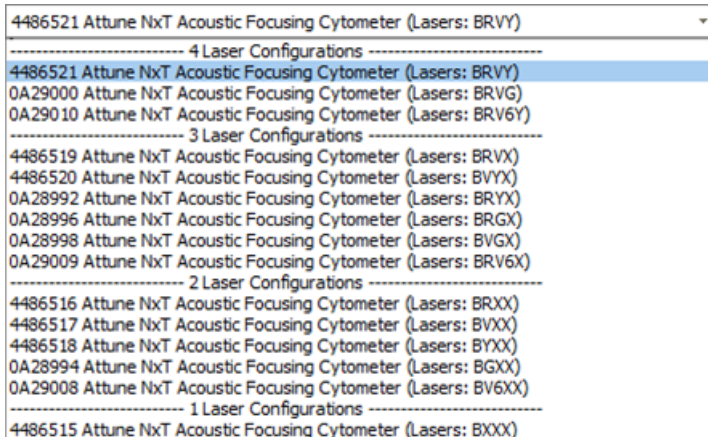
The screenshot shows the 'Options' dialog box with the 'Configuration' tab selected. The main area is titled 'Access resources for Attune™ Cytometer Hardware' and contains the following sections:

- Hardware / Virtual Laser Configuration:**
  - Instrument Type: Attune CytPix Flow Cytometer
  - Attune Model: 0A48663 Attune CytPix Flow Cytometer (Lasers: BRVY)
  - Autosampler Model: CytKick™ Autosampler
- Enable Bubble Sensor:**
  - Enabled
- Automation Communication Settings:**
  - TCP/IP Communication Port: Valid ports are 1025-49151 (Default is 8225)
- Thermo Fisher Cloud Device Registration:**
  - Device Name: [Empty text box]
  - Register Device button

At the bottom right of the dialog are 'OK' and 'Cancel' buttons.

## Hardware/Virtual laser configuration

*Hardware/Virtual Laser Configuration* dropdown allows you select a virtual instrument configuration. This aids in the creation of Experiments when not connected to a cytometer by maintaining correct instrument settings and channel mapping.



- Clicking the **Hardware/Virtual Laser Configuration** combo box control opens the instrument configuration dropdown list, which is populated from the Attune™ NxT instrument configurations enumerated in the database.
- The available configurations are described in “Default instrument configurations” (“Default instrument configurations” on page 527).



## Overview

### Baseline calculation and Performance testing

To ensure that the Attune™ Cytometer is running in good condition, the instrument has a quality control feature referred to as *Performance Tracking*, which has two separate but related parts, **Baseline Calculation (BL)** and **Performance Test (PT)**.

On brand-new instruments, brand-new beadlots, or instruments that have gone through major service repair, a **Baseline Calculation** is performed. This is an in-depth quality control test that measures and sets the baseline standards for the instrument performance using Attune™ Performance Tracking Beads.

For the **Baseline Calculation**, the instrument measures the median fluorescence intensity of each bead and the r%CV (robust percent coefficient of variation) in all fluorescence detectors to determine cytometer settings and provide target values for subsequent application-specific settings (see “Baseline calculation and Performance test data” on page 540). The **Baseline Calculation** also includes the measurement of **PMT Functional Response**, which measures and calculates the optimal voltages of the instrument per fluorescent parameter (see “Baseline Functional Response (BFR)” on page 540).

After the Baseline values are established, the same lot of Attune™ Performance Tracking Beads are used to run the daily **Performance Test**. The **Performance Test** determines variation from the Baseline values to track the daily performance of the cytometer, where deviations in performance can indicate a need for maintenance or servicing.

### Performance tracking workflow (when to run)

#### Baseline Calculation

Baseline Calculation is performed:

- By the Field Service Engineer (FSE) at time of installation.
- By the FSE after any major service.
- By the user every time the performance tracking bead lot changes.
- By the user when recommended by FSE or FAS.

#### Performance Test

Performance Test is performed by the user every day samples are run or recorded.

---

**Note:** **Baseline Calculation** can only be performed by Administrator or Advanced User accounts. **Performance Test** can be performed by Administrator, Advanced User, and User accounts.

---

## Baseline calculation and Performance test data

**Baseline calculation** uses Attune™ Performance Tracking Beads to generate data for the parameters listed below to establish the **initial** status of the cytometer. The results are reported in the **Baseline Test Results** and the **Baseline Functional Response** screens.

- PMTV (photomultiplier tube voltage) are adjusted to place the brightest bead at target MFI (median fluorescence intensity) values; voltage for each channel is reported.
- Target MFI (target median fluorescence intensity) is reported.
- Measured MFI is determined and reported.
- r%CV (robust percent coefficient of variation) of the brightest bead is recorded.
- Relative Quantum Efficiency (Qr) of each detector is recorded.
- Relative Background (Br) of each detector is determined.
- Linear regression (Linearity) is calculated and recorded.
- Area Scaling Factor (ASF) is calculated and reported for every laser and automatically updated in **Advanced Settings**.
- Laser Delay setting is automatically calculated.

**Performance test** uses Attune™ Performance Tracking Beads to determine the variation of the test data from the target values established by the Baseline calculations to track the daily performance of the instrument. The results are reported in the **Performance Test Results** screen.

- Same processes and measurements as Baseline are determined and reported.
- Delta PMTV (change in PMT voltage) from Baseline is also reported.

## Baseline Functional Response (BFR)

To obtain high quality fluorescent flow cytometer data, a well-optimized instrument is required. To identify optimal voltages for each detector, users must spend 1–2 days performing an experiment known as a "voltage walk" or "voltration". This requires calibration beads or multiple samples stained with conjugated antibody fluorophores to be recorded, gated, and analyzed to find the desired voltage for each of the 14 fluorescent detectors.

The **Baseline Functional Response (BFR)** function of the Attune™ Cytometric Software, performed automatically by the instrument before **Baseline Calculation**, determines the **Optimal PMTV** for each fluorescent detector in 10–20 minutes and records them in the **Baseline Test Results** (see "Completion of Baseline" on page 547). This enables users to optimize the instrument for their experiments by suggesting the minimum voltage required for the best signal more efficiently to ensure that:

1. Negative populations are above the noise on the low end of the measurement.
2. The positive populations are below the top of the linear range at the high end of the measurement.

---

**Note:** The optimal voltages are only calculated during the **Baseline** run. Running a **Performance Test** does not calculate optimal voltages. **Performance Tests** can be performed after a **Baseline** has been successfully performed. It is not necessary to perform extra **Baseline Calculations** outside of

the above listed recommendations due to the stability of the instrument, which can be monitored by observing the low delta PMTV values recorded during the **Performance Test**.

**IMPORTANT!** Use the **BFR** as a starting point for the optimization of the instrument. Voltages should be fine tuned for any application for brighter signals in the red channel.

## Baseline and Performance Test (BL/PT) module

The **Baseline and Performance Test (BL/PT) module** of the Attune™ Cytometric Software is used to:

- “Set up and run a Baseline” on page 542 (includes the BFR measurement)
- “Run a Performance test” on page 548
- Chapter 20, “Performance Test reports”

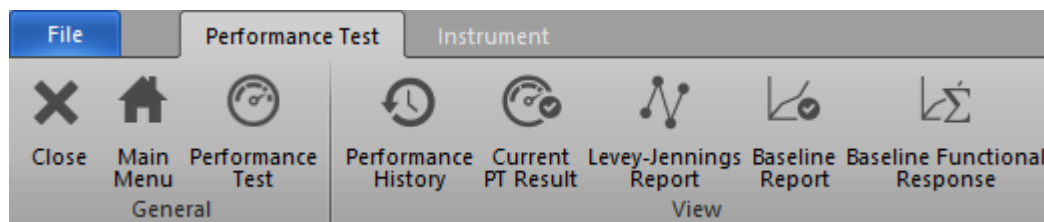
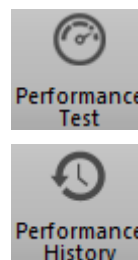
**Note:** For information about viewing the **Baseline** and **Performance Test reports**, see Chapter 20, “Performance Test reports”.

### Open the BL/PT module

On the **Main Menu** (“Main Menu” on page 46), click **Performance Test** to open the **BL/PT module**, which contains the **Performance Test** ribbon tab.

Alternatively, click **Performance History** on the **Instrument** ribbon tab (“Instrument tab” on page 88), when the instrument is connected.

- In the **BL/PT module**, only the **File**, **Home**, **Performance Test**, and **Instrument** ribbon tabs are available. All other tabs are hidden.



**Note:** The **Performance Test** ribbon tab is only visible when the **BL/PT module** is active.

- If no **Baseline** has previously been run, the **Baseline setup** screen is displayed; see “Set up and run a Baseline” on page 542.
- If a **Baseline** already exists, the **Performance Test setup** screen is displayed; see “Run a Performance test” on page 548.
- To close the **BL/PT module** and return to the previous view, click **Close**.
- To return to the **Main Menu**, click **Main Menu**.

# Set up and run a Baseline

## Baseline setup screen

Open the **BL/PT module** (“Baseline and Performance Test (BL/PT) module” on page 541) to view the **Baseline setup** screen.

---

**Note:** If a Baseline already exists, the **Performance Test setup** screen is displayed instead of the **Baseline setup** screen. On the **Performance Test setup** screen (“Run a Performance test” on page 548), you can run a Performance Test or reset the Baseline.

---

The **Baseline setup** screen provides general instructions for setting up a Baseline:

1. Check instrument configuration.
2. Verify bead lot number.
3. Add 4 drops bead solutions to 3 mL of Focusig Fluid.
4. Load tube.

For detailed procedures, see “Set up and run a Baseline” on page 542.

---

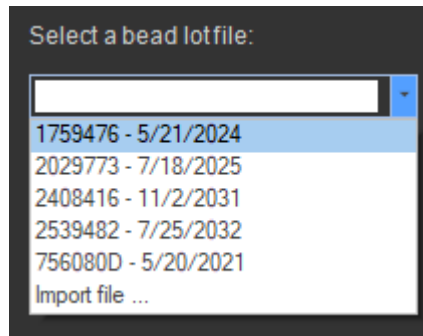
**Note:** **Baseline Calculation** can only be performed by users with Administrator or Advanced User accounts.

**Performance Test** can be performed by Administrator, Advanced User, and User accounts.

---



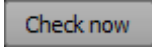
- The **Select bead lot file** dropdown list lists all bead lot files that have been imported.



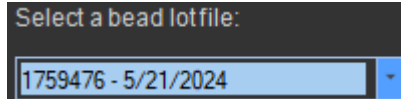
- Each file in the dropdown list is listed in the following format:  
**L/N: ###** = Bead lot number  
**DD/MM/YYYY** = Expiration date  
 The bead lot number is the **first seven digits** printed on the label (ignore the alpha numeric characters).
- To set up a **Baseline** for another bead lot, select the bead lot file of interest from the dropdown list. The software displays the bead lot number and expiration date for the currently selected bead lot.

## Set up and run a baseline

1. Click **Check now** to verify that the current filter configuration matches the default instrument configuration. If there is a mismatch between the current filters and the default instrument configuration, there is a high possibility of Baseline or Performance test failure.

A rectangular button with a light gray background and a dark gray border, containing the text "Check now" in a dark gray font.

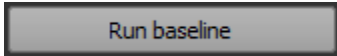
2. Select the bead lot file of interest from the **Select bead lot file** dropdown list.



3. Verify that the bead lot number you have selected agrees with the bead lot number of the Attune™ Performance Tracking Beads you are using.

The bead lot number is the **first seven digits** printed on the label (ignore the alpha numeric characters).

4. Prepare the beads by adding 4 drops of bead solution to 3 mL of Attune™ Focusing Fluid.
5. Load the tube as described in the *Attune™ NxT Flow Cytometer User Guide* (Part. No. [MAN0026547](#)) or the *Attune™ CytPix™ Flow Cytometer User Guide* (Pub. No. [MAN0019440](#)).
6. Click **Run Baseline**.

A rectangular button with a light gray background and a dark gray border, containing the text "Run baseline" in a dark gray font.

---

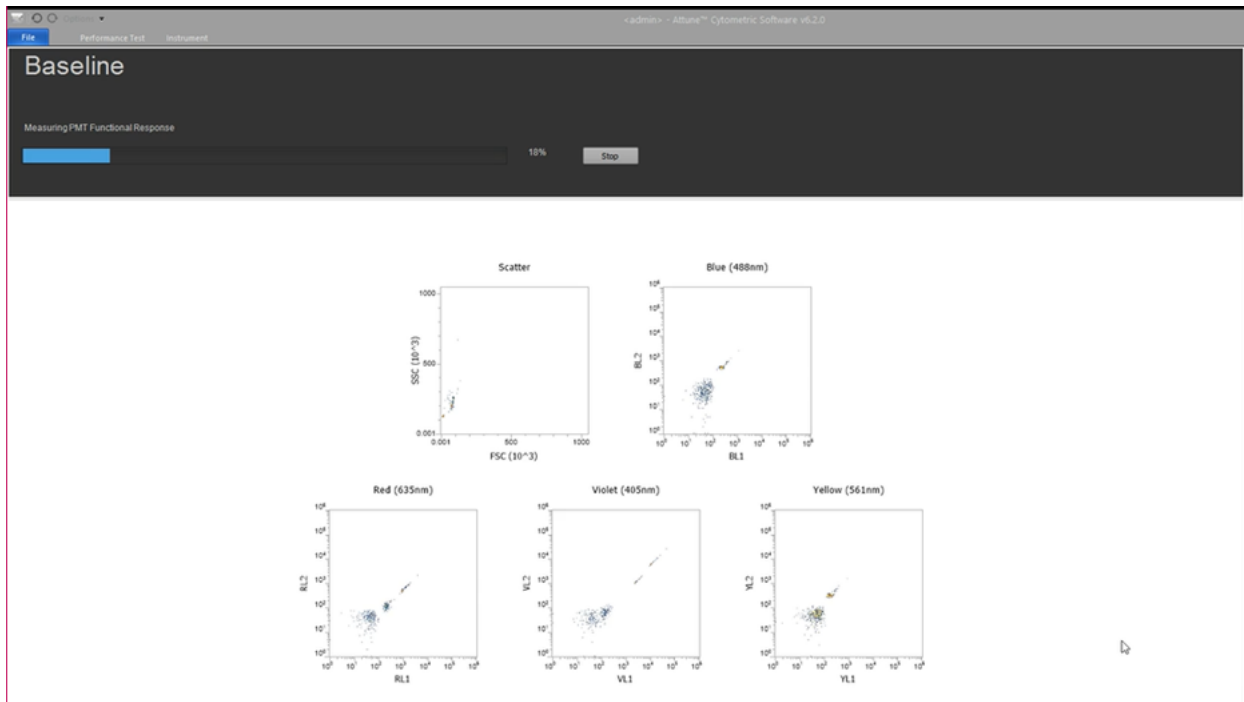
**Note:** If the **Startup** procedure has not been performed, the button displays **Run Startup** instead of **Run baseline**. You need to run the **Startup** procedure before proceeding.

---

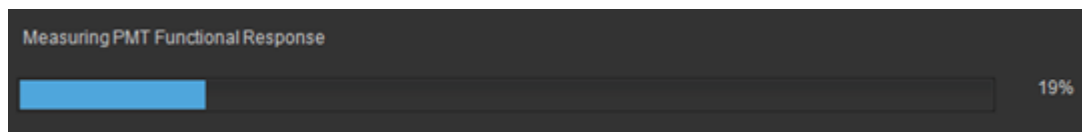
7. If the bead lot is expired, a warning dialog appears. Click **Yes** to continue or click **No** to cancel the Baseline run.  
If they are not already powered on, all lasers are powered on and remain on at the completion of the test process.

## Baseline screen

The **Baseline** screen provides progress information for the Baseline procedure.



- Five Baseline steps are displayed above the progress bar as they occur:
  - “Finding Beads”
  - “Measuring Laser Delays”
  - “Measuring PMT Functional Response”
  - “Adjusting PMTV to MFI Criteria”
  - “Gathering Final Statistics”
- Baseline procedure gathers statistics at flow rates of 100, 500, and 1000  $\mu\text{L}/\text{minute}$  and reports the results for 100  $\mu\text{L}/\text{minute}$  in the **Baseline Test Results** screen at the completion of the Baseline procedure.
- In the final step of the Baseline procedure, the instrument performs a fluid line calibration step.
  - Fluid line calibration requires a rinse to clear the fluid lines of sample.
  - Calibration compensates for fluid line volume tolerances and minimizes dead volume.
- The progress bar displays the percent completion (0–100%).



**Note:** The Baseline steps and the progress bar are updated by the software in real time. The BFR procedure corresponds to the 15–73% part of the Baseline progress.

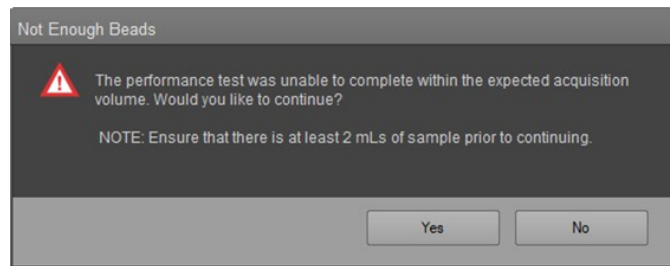
- Depending on the number of configured channels, the screen displays:
  - One scatter density plot (SSC-H vs. FSC-H)
  - Bivariate dot plots for each laser (for example, BL2-H vs. BL1-H)

For more information about the plots, see “Plots” (“Plots” on page 138).

- To zoom on a plot, use the **Size slider** on the **Application status bar** (“Size slider” on page 70). You cannot resize objects on the workspace. The plots are not resized when the application window is resized.
- To stop the Baseline run, click **Stop**. The baseline procedure stops and the software returns to **Baseline setup** (“Set up and run a Baseline” on page 542).



- If the sample is exhausted before the Baseline completes, the software displays the **Not Enough Beads** dialog, which prompts the user to add more solution to enable the instrument to draw more sample to complete the Baseline.



- At the completion of the Baseline procedure, the **Baseline Test Results** screen is displayed (see “Completion of Baseline” on page 547).



# Completion of Baseline

## Baseline completion without errors

If the Baseline completes without errors, the software displays the **Baseline Test Results** with the **Update Default Instrument Settings** dialog.

The screenshot shows the 'Baseline Test Results' screen. At the top, it indicates 'Baseline test successful' with a green checkmark. Below this, there is a table with columns: Channel, PMTV, Optimal PMTV, Target MFI, MFI, Robust %CV, Qr, Background, Linearity, ASF, Laser Delay, and Result. The table lists various channels (FSC, SSC, BL1-3, RL1-3, VL1-4, YL1-4) with their respective values. A dialog box titled 'Update Default Instrument Settings' is overlaid on the table, asking 'Optimal PMTV values have been calculated on the system. Would you like to update your default voltage settings with calculated values?' with 'Yes' and 'No' buttons.

Channel	PMTV	Optimal PMTV	Target MFI	MFI	Robust %CV	Qr	Background	Linearity	ASF	Laser Delay	Result
FSC	369	--	300,000	298,705	2.24 %	0.000	0	0.000	1.00	1100	✓
SSC	275	--	300,000	301,083	4.42 %	0.000	0	0.000	1.00	1100	✓
BL1	420	403	300,000	295,946	1.63 %	0.037					✓
BL2	378	394	300,000	295,983	1.50 %	0.042					✓
BL3	457	461	300,000	300,682	2.97 %	0.015					✓
RL1	355	367	300,000	297,420	2.98 %	0.076					✓
RL2	381	382	300,000	300,970	2.97 %	0.011					✓
RL3	434	410	300,000	297,070	3.22 %	0.046					✓
VL1	249	289	300,000	300,818	0.83 %	0.019	3067	1.000	1.10	701	✓
VL2	291	293	300,000	295,296	1.01 %	0.015	1033	0.997	1.10	701	✓
VL3	373	416	300,000	297,070	1.48 %	0.018	93	1.000	1.10	701	✓
VL4	398	414	300,000	299,522	2.65 %	0.002	283	0.990	1.10	701	✓
YL1	436	438	300,000	297,844	1.52 %	0.109	59	1.000	0.91	313	✓
YL2	410	423	300,000	296,356	2.50 %	0.068	38	1.000	0.91	313	✓
YL3	468	456	300,000	296,884	3.09 %	0.013	197	1.000	0.91	313	✓
YL4	475	479	300,000	297,090	3.58 %	0.005	369	1.000	0.91	313	✓

- The **Baseline Test Results** represents initial status of the cytometer and displays the data described in “Baseline calculation and Performance test data” on page 540, including Optimal PMTV values. For more information about the **Baseline Test Results** and how to interpret them, see “Baseline Report” on page 572.
- The **Update Default Instrument Settings** dialog asks if the user would like to update or overwrite their default system instrument voltages for the optimal voltages calculated in the Baseline.
  - Click **Yes** to use the optimal settings calculated in the Baseline in all newly created Experiments. This does not change templates or existing Experiments.
  - Click **No** if you do not want to use the optimal values in newly created Experiments.

## Baseline completion with errors

If the Baseline completes with errors, the software displays the **Errors Detected** dialog, which provides a list of the channel statistics that failed.

To go to the **Baseline Test Results** screen (“Baseline Report” on page 572), click **Continue to results**.

## Run a Performance test

### Performance test setup screen

Open the **BL/PT module** (“Baseline and Performance Test (BL/PT) module” on page 541) to view the **Performance Test setup** screen.

---

**Note:** If no Baseline exists, the **Baseline setup screen** is displayed instead of the **Performance Test setup** screen. You need to run the Baseline first; see “Set up and run a Baseline” on page 542.

---

The **Performance Test setup** screen provides general instructions for setting up a Performance Test:

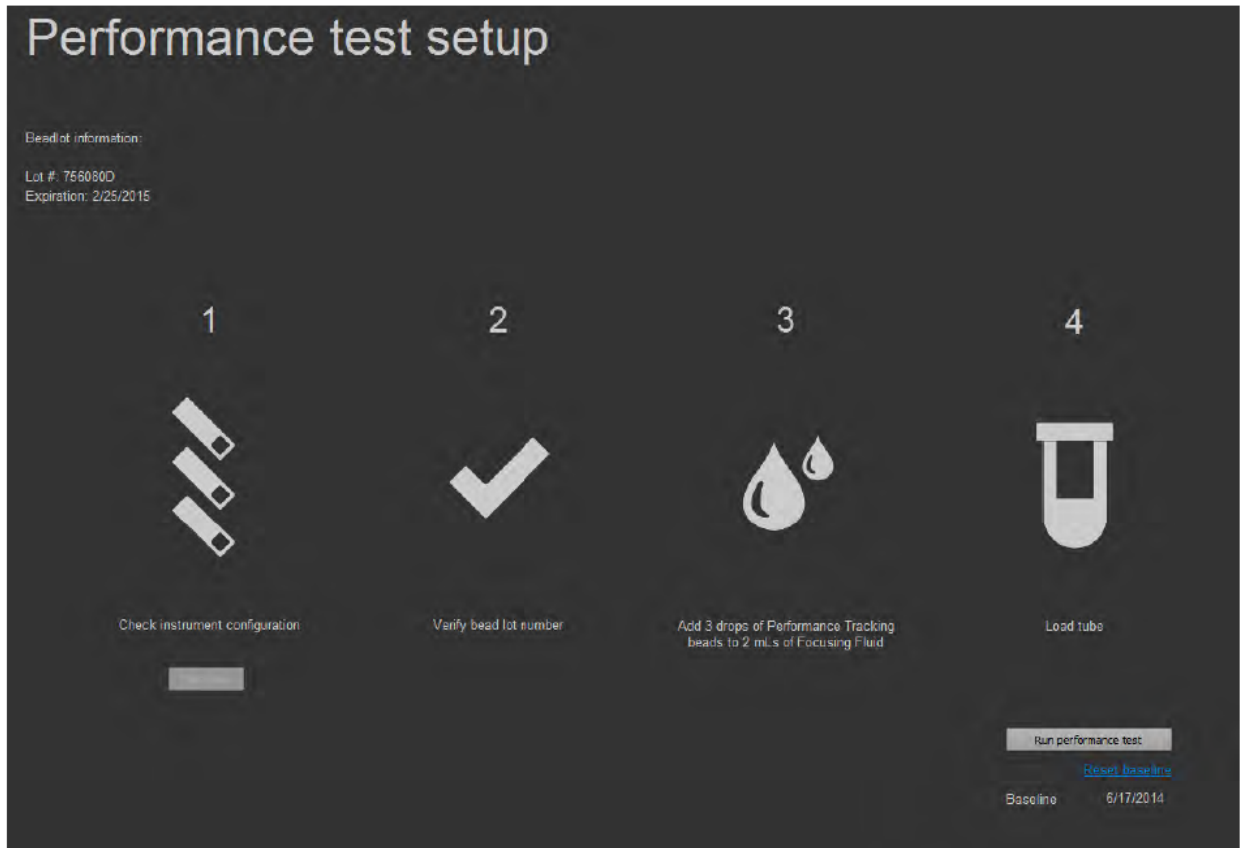
1. Check instrument configuration.
2. Verify bead lot number.
3. Add 3 drops bead solution to 2 mL Focusing Fluid.
4. Load tube.

For detailed procedures, see “Set up and run a Performance test” on page 550.

---

**Note:** **Baseline Calculation** can only be performed by users with Administrator or Advanced User accounts.

**Performance Test** can be performed by Administrator, Advanced User, and User accounts.



- The software displays the bead lot number and expiration date for the most recent bead lot used to set up a Baseline in the following format:

Lot #: 2029773  
Expiration: 12/30/2023

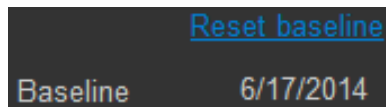
**L/N: ###** = Bead lot number

**DD/MM/YYYY** = Expiration date

The bead lot number is the **first seven digits** printed on the label (ignore the alpha numeric characters).

- If the bead lot has expired, the lot number and expiration date are in contrasting color.

- The software displays the current Baseline information in the lower right of the screen. To reset the current Baseline, click **Reset Baseline**, which opens the **Baseline setup** screen (“Set up and run a Baseline” on page 542).



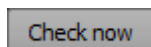
---

**Note:** The **Reset Baseline** hyperlink is active only if you are assigned the privileges to run the Baseline (see “Account permissions” on page 44).

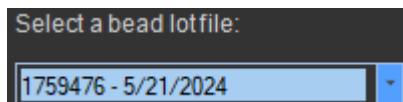
---

## Set up and run a Performance test

1. Click **Check now** to verify that the current filter configuration matches the default instrument configuration. If there is a mismatch between the current filters and the default instrument configuration, there is a high possibility of Baseline or Performance test failure.



2. Select the appropriate bead lot file from the **Select bead lot file** dropdown list. For detailed procedures, see “Set up and run a Baseline” on page 542.



If needed, you can first select a different Baseline to run the Performance Test against, then click **Reset Baseline** to open the **Baseline setup** screen.

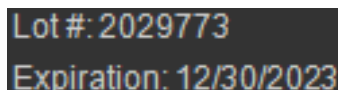


---

**Note:** The **Reset Baseline** hyperlink is active only if you are assigned the privileges to run the Baseline (see “Account permissions” on page 44).

---

3. Verify that the bead lot number you have selected agrees with the bead lot number of the Attune™ Performance Tracking Beads you are using.



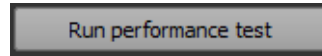
The bead lot number is the **first seven digits** printed on the label (ignore the alpha numeric characters).

4. Prepare the test beads by adding 3 drops of bead solution to 2 mL of Attune™ Focusing Fluid.

5. Load the tube to the cytometer.

For more information, see the *Attune™ NxT Flow Cytometer User Guide* (Part. No. [MAN0026547](#)) or the *Attune™ CytPix™ Flow Cytometer User Guide* (Pub. No. [MAN0019440](#)).

6. Click **Run Performance Test**.



---

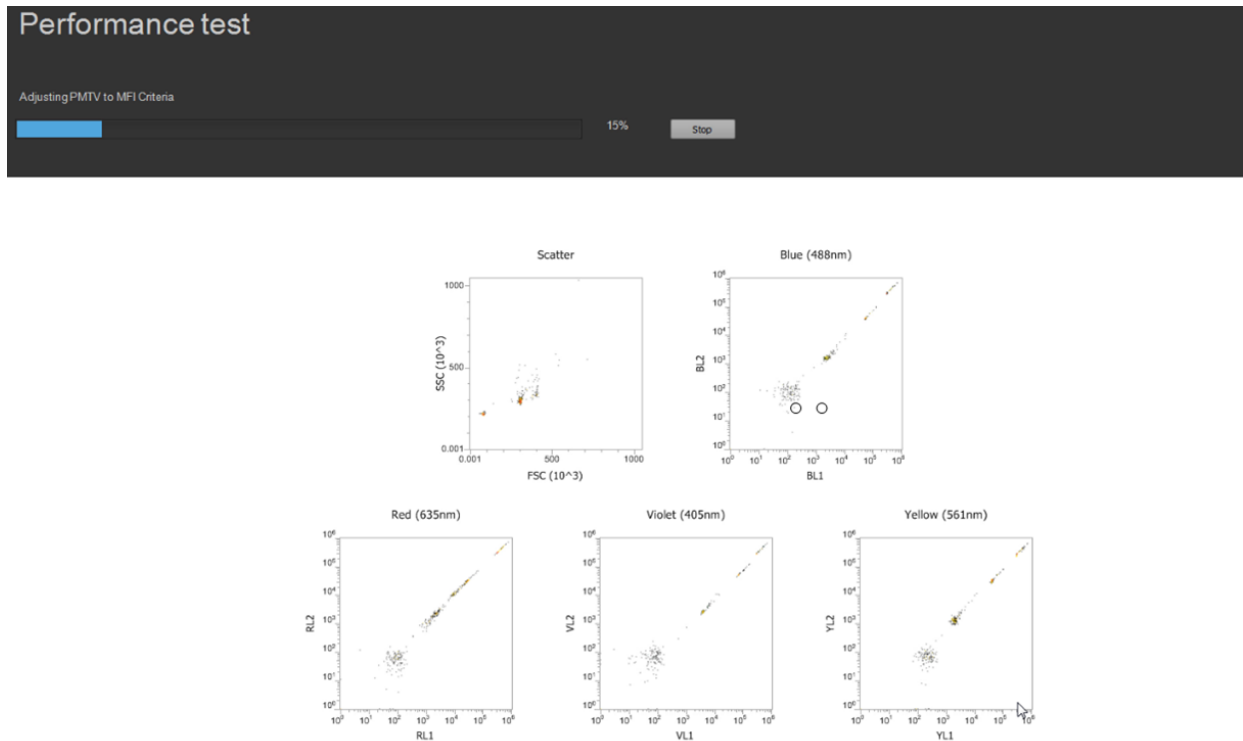
**Note:** If the **Startup** procedure has not been performed, the button displays **Run Startup** instead of **Run Performance Test**. You need to run the **Startup** procedure before proceeding.

---

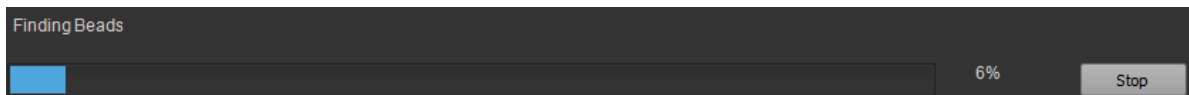
7. If the bead lot is expired, a warning dialog appears. Click **Yes** to continue or click **No** to cancel the Performance Test run.  
If they are not already powered on, all lasers are powered on and remain on at the completion of the test process.
8. After the Performance Test is complete, click **Sanitize Attune SIP** (“Functions group” on page 90) to run the **SIP sanitize** procedure before proceeding with the experiments.

## Performance test screen

The **Performance test** screen provides progress information for the Performance Test procedure.



- Four Performance Test steps are displayed above the progress bar as they occur:
  - “Finding Beads”
  - “Measuring Laser Delays”
  - “Adjusting PMTV to MFI Criteria”
  - “Gathering Final Statistics”
- The progress bar displays the percent completion (0–100%).



**Note:** The Performance Test steps and progress bar are updated by the Attune™ Cytometric Software in real time.

- Depending on the number of configured channels, the screen displays:
  - One scatter density plot (SSC-H vs. FSC-H)
  - Bivariate dot plots for each laser (for example, BL2-H vs. BL1-H)

For more information about the plots, see “Plots” (“Plots” on page 138).

- To zoom on a plot, use the **Size slider** on the **Application status bar** (“Size slider” on page 70). You cannot resize objects on the workspace. The plots are not resized if the application window is resized.
- To stop the Performance Test procedure, click **Stop**. The procedure stops and the software returns to Baseline setup (“Set up and run a Baseline” on page 542).



## Completion of Performance test

### Performance test completion without errors

If the Performance Test completes without errors, the software:

- Displays the **Performance Test Results** screen (“Current PT results” on page 561).

**Performance Test Results**

✓ Performance test successful

Baseline: 1759476 - 4/3/2017

6/8/2017 6:41:55 PM

Channel	PMTV	Delta PMTV	Target MFI	MFI	Robust %CV	Qr	Background	Linearity	ASF	Laser Delay	Result
FSC	395	-4	300,000	302,628	2.23 %	0.000	0	0.000	1.06	1100	✓
SSC	360	0	300,000	290,538	2.94 %	0.000	0	0.000	1.06	1100	✓
BL1	417	-2	300,000	301,217	1.49 %	0.060	132	1.000	1.06	1100	✓
BL2	352	-3	300,000	305,144	1.17 %	0.054	177	1.000	1.06	1100	✓
BL3	437	-6	300,000	301,522	2.44 %	0.039	22	0.999	1.06	1100	✓
RL1	380	-4	300,000	300,287	3.58 %	0.003	13	0.981	1.03	1415	✓
RL2	371	-3	300,000	306,141	3.54 %	0.000	43	0.947	1.03	1415	✓
RL3	392	-3	300,000	304,672	3.57 %	0.004	23	0.948	1.03	1415	✓
VL1	322	1	300,000	296,147	1.35 %	0.018	1350	0.999	1.08	655	✓
VL2	374	1	300,000	296,768	1.66 %	0.022	419	0.996	1.08	655	✓
VL3	374	-1	300,000	302,433	2.57 %	0.030	76	1.000	1.08	655	✓
VL4	423	-3	300,000	300,499	3.19 %	0.004	264	0.999	1.08	655	✓
YL1	375	-2	300,000	306,149	1.55 %	0.117	94	0.999	1.01	329	✓
YL2	402	0	300,000	294,123	2.81 %	0.072	27	1.000	1.01	329	✓
YL3	406	-2	300,000	296,638	4.00 %	0.008	147	0.996	1.01	329	✓
YL4	467	-2	300,000	299,562	4.51 %	0.003	306	0.993	1.01	329	✓

- Calculates the system area scaling factor constant based on the Performance Test results for each configured laser.
- Updates the default system values for the area scaling factor in the Instrument Settings panel (“Area scaling factor (ASF)” on page 400).
- Updates the laser delay.

---

**IMPORTANT!** The optimal voltages are only calculated during the Baseline Functional Response (BFR) part of the Baseline procedure. Performance Test does not calculate optimal voltages.

Voltage settings are not updated or modified as a result of daily Performance testing. **For best results, use BFR settings as starting points and manually fine tune voltage setting for each experiment.** Laser delays for each flow rate are set based on the Performance Test results. These values cannot be modified.

---

## Performance test completion with errors

If the Performance Test completes with errors, the software displays the **Errors Detected** dialog, which provides a list of the channel statistics that failed

To go to the **Performance Test Results** screen (“Current PT results” on page 561), click **Continue to results**.

## Performance test reports

After running the Performance Tests, you can generate the following reports in the Attune™ Cytometric Software:

- Performance History
- Current PT Results
- Levey-Jennings Report
- Baseline Report
- Baseline Functional Response

For information about viewing the Performance Test reports, see Chapter 20, “Performance Test reports”.







## Overview

**Performance test** is a comprehensive set of procedures to monitor the daily performance of the cytometer. The **Performance test status** for the day is displayed below the Sign in text fields on the **Login screen**.

The **Performance test status** includes a message and an icon that indicate the date and the results of the last performance test (see below). If no baseline has ever been run, no message is displayed.

**Note:** **Performance test status** is also displayed on the **Main Menu** below the Main Menu buttons.

Icon	Login Screen	Main Menu
	Performance test has been completed successfully today.	Performance test has been completed successfully today.
	Performance test has not been completed today.	Performance test has not been completed today.
	Performance test has not been completed successfully today.	Performance test has not been completed successfully today.
		Bead lot has expired (see <b>Note</b> below) <sup>[1]</sup>

<sup>[1]</sup> If the baseline calculations have expired, the performance test status bar displays the “Bead lot has expired” message. For more information on baseline calculations, see “Set up and run a Baseline” on page 542, and “Completion of Baseline” on page 547.

## Performance tests

The Attune™ Cytometric Software uses Attune™ Performance Tracking Beads to define the cytometer’s **Baseline status**. During this process, the median fluorescence intensity of each bead and the r%CV (robust percent coefficient of variation) are measured in all fluorescence detectors. Software algorithms use this information to determine cytometer settings and provide target values for subsequent application specific settings (see “Performance Test data”, below).

After the **Baseline** values are established, you use the same lot of Attune™ Performance Tracking Beads to run the **Performance Tests**. The **Performance Tests** measure variation from the **Baseline** values to track the daily performance of the cytometer.

**Note:** For information about running **Performance Tests**, see Chapter 19, “Performance testing”.

## Baseline calculation and Performance test data

**Baseline calculation** uses Attune™ Performance Tracking Beads to generate data for the parameters listed below to establish the **initial** status of the cytometer. The results are reported in the **Baseline Test Results** and the **Baseline Functional Response** screens.

- PMTV (photomultiplier tube voltage) are adjusted to place the brightest bead at target MFI (median fluorescence intensity) values; voltage for each channel is reported.
- Target MFI (target median fluorescence intensity) is reported.
- Measured MFI is determined and reported.
- r%CV (robust percent coefficient of variation) of the brightest bead is recorded.
- Relative Quantum Efficiency (Qr) of each detector is recorded.
- Relative Background (Br) of each detector is determined.
- Linear regression (Linearity) is calculated and recorded.
- Area Scaling Factor (ASF) is calculated and reported for every laser and automatically updated in **Advanced Settings**.
- Laser Delay setting is automatically calculated.

**Performance test** uses Attune™ Performance Tracking Beads to determine the variation of the test data from the target values established by the Baseline calculations to track the daily performance of the instrument. The results are reported in the **Performance Test Results** screen.

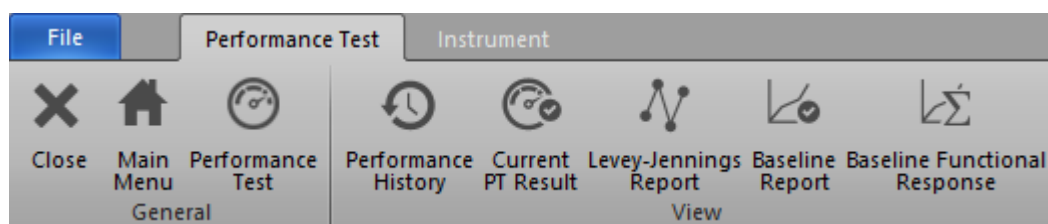
- Same processes and measurements as Baseline are determined and reported.
- Delta PMTV (change in PMT voltage) from Baseline is also reported.

## Performance test reports

After running the **Performance Tests**, you can generate the following reports in the Attune™ Cytometric Software:

- Performance History Report (“Performance History Report” on page 557)
- Current PT Results (“Current PT results” on page 561)
- Levey-Jennings Report (“Levey-Jennings Report” on page 565)
- Baseline Report (“Baseline Report” on page 572)
- Baseline Functional Response (“Baseline Functional Response (BFR) Report” on page 576)

To view a report, click the corresponding button on the **Performance Test** tab of the **BL/PT module**.

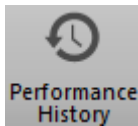


## Performance History Report

The **Performance History Report** shows the pass/fail status of all **Performance Tests** run against a selected **Baseline test** during a selected time period. The **Performance History Report** is useful for gauging the overall “health” of the instrument over the selected time period.

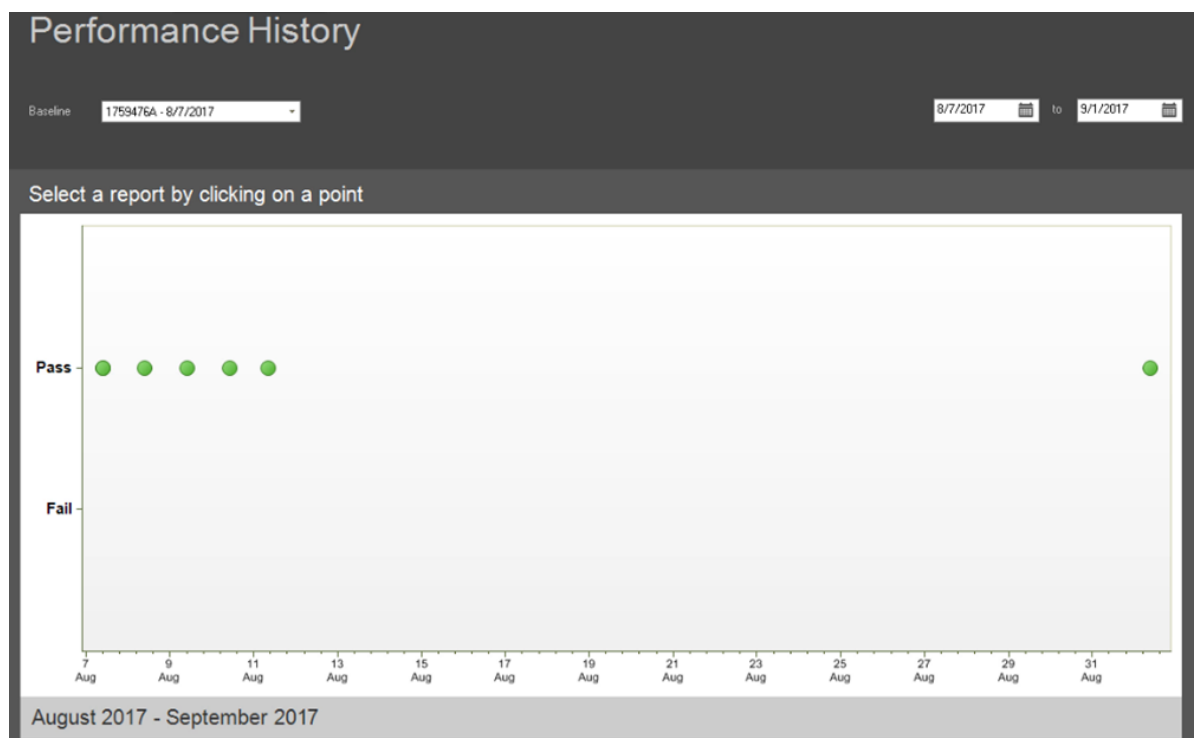
### View the Performance History Report

On the **Performance Test** ribbon (available on the **BL/PT module**; “Baseline and Performance Test (BL/PT) module” on page 541), click **Performance History**.

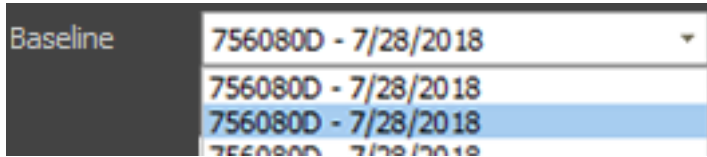


Alternatively, click **Performance History** on the **Instrument** ribbon tab (“Instrument tab” on page 88).

- By default, the **Performance History Report** is displayed for the most recent **Baseline**, which shows the **Pass/Fail status** of **Performance Tests** for the most recent 180-day period. For more information, see “Results plot” on page 559.



- To view the report for **Performance Tests** associated with another Baseline and Bead lot, select the Baseline of interest from the **Baseline** dropdown list.



Each Baseline in the dropdown list is listed in the following format:

**###** = Performance Test bead lot number

**DD/MM/YYYY** = date Baseline status was established

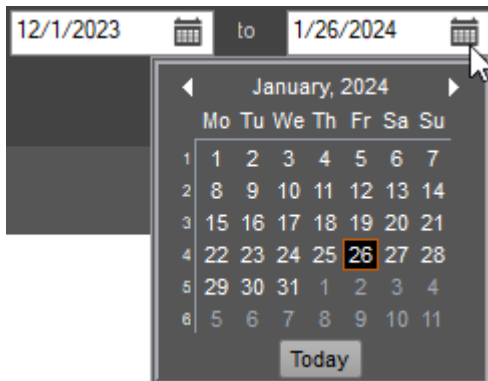
---

**Note:** The **Baseline** dropdown list is enabled only if more than one Baseline test has been run. The Baselines are ordered chronologically, with the most recent displayed first.

---

- By default, the **Performance History Report** shows the most recent 6-month data for the **Performance Test** series in the selected **Baseline**.

To view the **Performance History Report** for a different data range, click the **calendar** icon on the **Date Range** tool to select the desired beginning and end dates.




---

**Note:** By default, the range of the **Performance History Report** begins with the first **Baseline** and ends with the most recent **Performance Test** available for the selected **Baseline**.

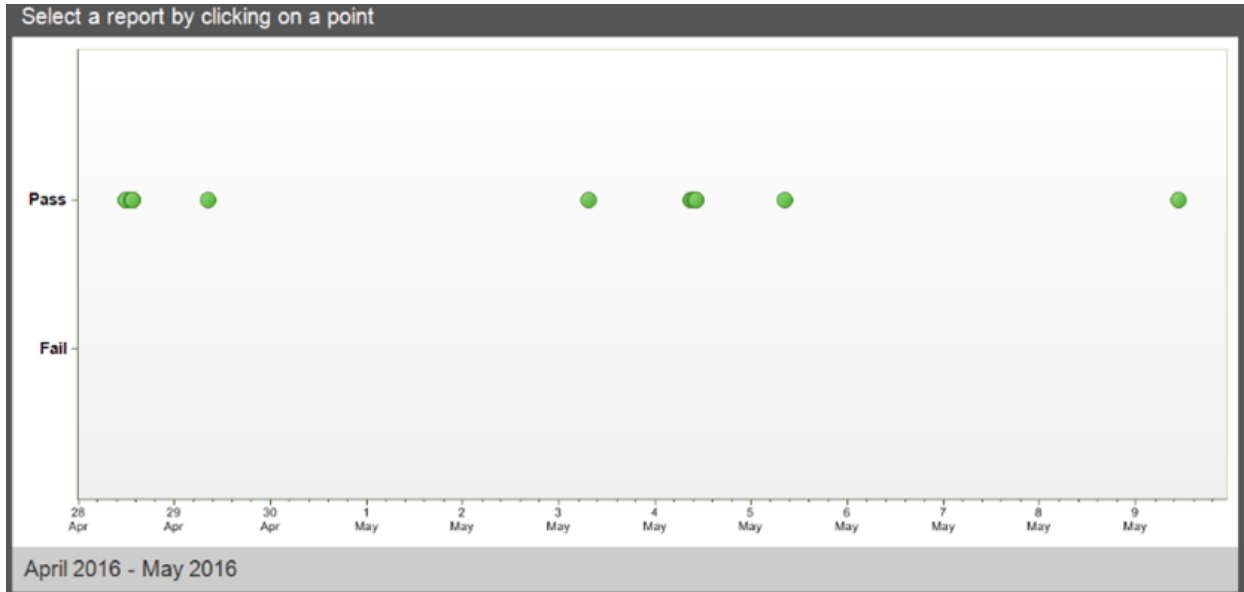
If the default range exceeds 180 days, then the report shows data for the most recent 180-day time period.

Using the **Date Range** tool, you can select a date range of up to 180 days.



---

## Results plot

The **Results plot** graphically summarizes the Pass/Fail status of the selected Baseline and associated Performance Tests within the selected date range.



- Each data point on the **Results plot** corresponds to a single date and the color of the data point indicates whether the result was a pass or fail.

Icon	Pass/Fail status
	<b>Pass:</b> All Performance Tests run on that date are within the limits set by the Baseline calculation.
	<b>Fail:</b> One or more Performance Tests run on that date deviate from the target set by the Baseline calculation.

- If more than one result is available on a single date and the range is set such that the results overlap, the most recent result is displayed first.
- Place the pointer over a data point to show the **Pass/Fail status** and the date of the Performance Test.
- Click a data point in the plot to open the Baseline or Performance Test result for that date.
- Use the mouse wheel to zoom in on the X-axis, which displays a scroll bar when the zoom function is applied. The Y-axis cannot be zoomed.
- If the date range is changed, the zoom is reset to the default.

**Note:** The **Pass/Fail status** is based on the overall result of the Performance Tests for that date and the Baseline.

- If all Performance Tests run on that date meet the target values as determined by the Baseline test, the status is **Pass**.
- If one or more Performance Tests run on that date do not meet the Baseline test values, the status is **Fail**.

## Print the performance history report

Select **File ▶ Print** to print the Performance History Report.

In addition to the Results plot, the printout includes a header containing the date range of the report and the instrument information.

Attune™ NxT Cytometer Baseline Report

7/28/2018 8:28:00 PM

---

Cytometer: Attune NxT Acoustic Focusing Cytometer (Lasers: BRVY) User: admin	
Instrument Model Number: 4486521	Institution: Thermo Fisher Scientific
Instrument Serial Number: 1AFC200010913	Software: 2.7.873.0

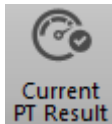
## Current PT results

**Current PT Results** shows the results of the current Performance Test and includes all statistics and calculations for all configured channels on the instrument.

### View the Current PT Results

**Current PT Results** report appears automatically after a Performance Test is completed.

You can also access the most recent Performance Test report by clicking **Current PT Result** button on the **Performance Test** ribbon tab of the **BL/PT module** (“Baseline and Performance Test (BL/PT) module” on page 541).



**Performance Test Results**



✓ Performance test successful

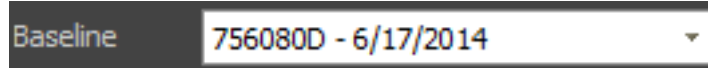
Baseline: 1759476 - 4/3/2017

6/8/2017 6:41:55 PM

Channel	PMTV	Delta PMTV	Target MFI	MFI	Robust %CV	Qr	Background	Linearity	ASF	Laser Delay	Result
FSC	395	-4	300,000	302,628	2.23 %	0.000	0	0.000	1.06	1100	✓
SSC	360	0	300,000	290,538	2.94 %	0.000	0	0.000	1.06	1100	✓
BL1	417	-2	300,000	301,217	1.49 %	0.060	132	1.000	1.06	1100	✓
BL2	352	-3	300,000	305,144	1.17 %	0.054	177	1.000	1.06	1100	✓
BL3	437	-6	300,000	301,522	2.44 %	0.039	22	0.999	1.06	1100	✓
RL1	380	-4	300,000	300,287	3.58 %	0.003	13	0.981	1.03	1415	✓
RL2	371	-3	300,000	306,141	3.54 %	0.000	43	0.947	1.03	1415	✓
RL3	392	-3	300,000	304,672	3.57 %	0.004	23	0.948	1.03	1415	✓
VL1	322	1	300,000	296,147	1.35 %	0.018	1350	0.999	1.08	655	✓
VL2	374	1	300,000	296,768	1.66 %	0.022	419	0.996	1.08	655	✓
VL3	374	-1	300,000	302,433	2.57 %	0.030	76	1.000	1.08	655	✓
VL4	423	-3	300,000	300,499	3.19 %	0.004	264	0.999	1.08	655	✓
YL1	375	-2	300,000	306,149	1.55 %	0.117	94	0.999	1.01	329	✓
YL2	402	0	300,000	294,123	2.81 %	0.072	27	1.000	1.01	329	✓
YL3	406	-2	300,000	296,638	4.00 %	0.008	147	0.996	1.01	329	✓
YL4	467	-2	300,000	299,562	4.51 %	0.003	306	0.993	1.01	329	✓

## Current PT Result report

- The overall **Pass/Fail status** of the current Performance Test is displayed by the **Pass/Fail indicator**, which includes a message and an indicator icon.  
**Pass:**  Performance test successful All channels meet the criteria set by the Baseline calculation.  
**Fail:**  Performance test completed with errors One or more of the channels fail to meet the criteria set by the Baseline calculation.
- To view the report for the Performance Tests associated with another Baseline and Bead lot, select the Baseline of interest from the **Baseline** dropdown list.

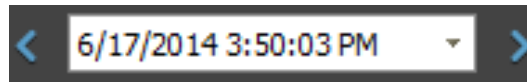


Each **Baseline** in the dropdown list is listed by the Performance Test bead lot number followed by the date the Baseline status was established

---

**Note:** The **Baseline** dropdown list is active only if more than one Baseline has been established. The Baselines are ordered chronologically, with the most recent displayed first.

- To view results from other Performance tests, select the desired test date from the **Performance Test history** dropdown, or click the blue **arrows** to navigate through the Baseline reports and Performance Test reports.



---

**Note:** If there are multiple reports for a single day, the reports are displayed chronologically, with the most recent displayed first.

---



## PT Results table


The **PT Results table** shows all statistics and calculations for all configured channels on the instrument.

Channel	PMTV	Target MFI	MFI	Robust %CV	Qr	Background	Linearity	ASF	Laser Delay	Result
FSC	381	300,000	298,275	1.90 %	0.000	0	0.000	1.04	1100	✓
SSC	316	300,000	293,378	3.84 %	0.000	0	0.000	1.04	1100	✓
BL1	467	300,000	298,406	1.19 %	0.049	112	1.000	1.04	1100	✓
BL2	417	300,000	297,482	1.05 %	0.045	158	1.000	1.04	1100	✓
BL3	473	300,000	294,423	1.81 %	0.033	40	1.000	1.04	1100	✓
RL1	393	300,000	298,334	3.62 %	0.115	60	0.999	1.00	1465	✓
RL2	421	300,000	292,456	3.60 %	0.012	328	0.994	1.00	1465	✓
RL3	478	300,000	298,215	3.94 %	0.058	152	0.997	1.00	1465	✓
VL1	286	300,000	296,710	1.42 %	0.016	1375	1.000	1.14	687	✓
VL2	313	300,000	298,217	1.70 %	0.021	552	0.998	1.14	687	✓
VL3	416	300,000	299,315	1.92 %	0.034	101	1.000	1.14	687	✓
VL4	437	300,000	297,411	2.55 %	0.004	270	0.993	1.14	687	✓
YL1	465	300,000	297,980	1.35 %	0.103	54	1.000	0.94	329	✓
YL2	446	300,000	299,817	1.28 %	0.083	30	1.000	0.94	329	✓
YL3	505	300,000	300,372	2.15 %	0.014	172	1.000	0.94	329	✓
YL4	514	300,000	295,153	2.87 %	0.006	301	1.000	0.94	329	✓

- The configured channels are listed in the following order and color-coded based on the laser color.

 : Scatter channels (FSC and SSC)

 : Blue channels

 : Red channels

 : Violet channels

 : Yellow channels

 : Green channels

- The data for the following parameters are displayed for all configured channels in the order specified below.
  - PMTV (photomultiplier tube voltage)
  - Delta PMTV
  - Target MFI (target median fluorescence intensity)
  - Measured MFI
  - Robust %CV (robust percent coefficient of variation)
  - Quantum Efficiency (Qr)
  - Background
  - Linearity
  - Area Scaling Factor (ASF)
  - Laser Delay

- The decimal places are only shown for **Robust %CV**, **Qr**, **Linearity**, and **Area Scaling Factor** (2 decimal places for Robust %CV and Area Scaling Factor, and 3 decimal places for Qr and Linearity).
- The **Pass/Fail status** for each configured channel is indicated by the status indicator icons shown below.
  - ✔ **Pass:** All statistics and calculations for the channel meet the criteria set by the Baseline calculation.
  - ⚠ **Fail:** One or more of the statistics or calculations for the channel deviate from the target set by the Baseline calculation.

---

**IMPORTANT!** The optimal voltages are only calculated during the **Baseline Functional Response (BFR)** part of the **Baseline** procedure. Running a **Performance Test** does not calculate optimal voltages, but still needs to be run daily. **Performance Tests** can be performed after a **Baseline** has been successfully performed. It is not necessary to perform extra **Baseline Calculations** outside of the recommended times due to the stability of the instrument (see “Performance tracking workflow (when to run)” on page 539), which can be monitored by observing the low delta PMTV values recorded during the **Performance Test**.

---

## Print the PT results report

Select **File** ▶ **Print** to print the PT Results Report.

In addition to the PT Results table, the printout includes a header containing the date range of the report, as well as the instrument information.

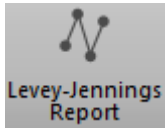
Cytometer: Attune NxT Acoustic Focusing Cytometer (Lasers: BRVG) User: admin	
Instrument Model Number: 0A29000	Institution: Thermo Fisher Scientific
Instrument Serial Number: 1AFC200010913	Software: 2.7.873.0

## Levey-Jennings Report

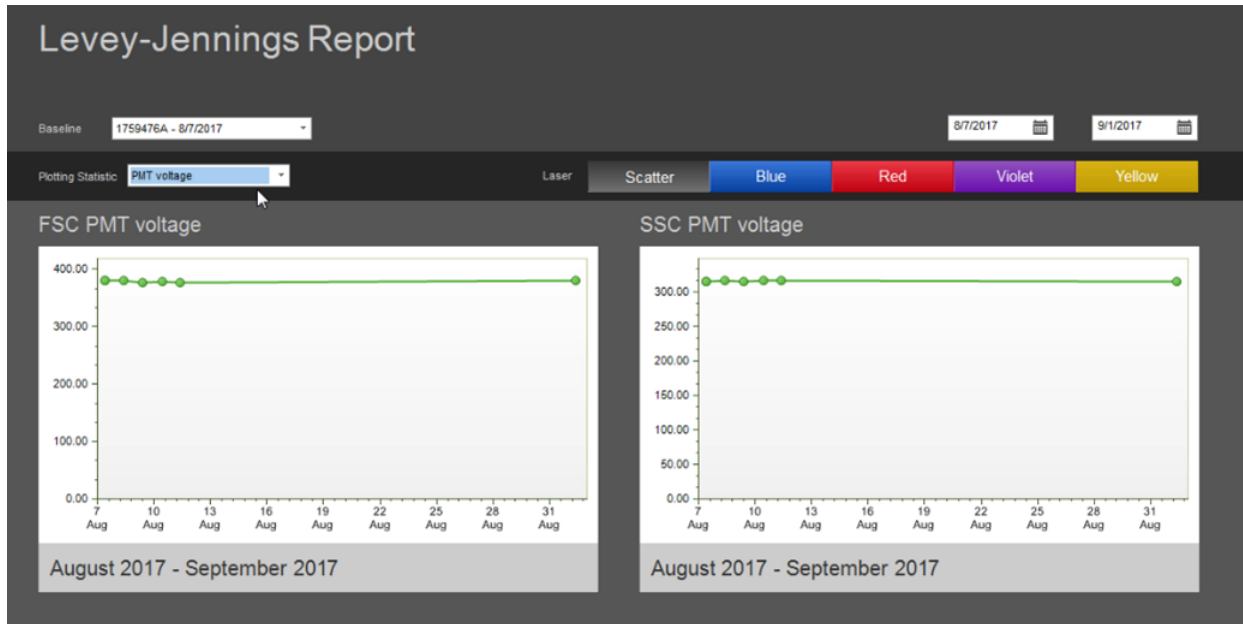
The **Levey-Jennings Report** tracks the parameters listed in “Baseline calculation and Performance test data” on page 556, for all configured channels to check for shifts and trends in cytometer performance and provides a visual indication of the cytometer performance over time.

### View the Levey-Jennings Report

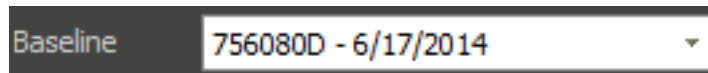
To view the Levey-Jennings Report, click **Levey-Jennings Report** on the **Performance Test** tab of the **BL/PT module** (“Baseline and Performance Test (BL/PT) module” on page 541).



The **Levey-Jennings Report** is shown for the most recent Baseline with data from the most recent 180-day period.



- To view the **Levey-Jennings Report** for Performance Tests associated with another Baseline and Bead lot, select the Baseline of interest from the **Baseline** dropdown list.



Each **Baseline** in the dropdown list is listed in the following format:

**###** = Performance Test bead lot number

**DD/MM/YYYY** = The date Baseline status was established

**Note:** The **Baseline** dropdown list is active only if more than one Baseline has been established. The Baselines are ordered chronologically, with the most recent displayed first.

- To view the **Levey-Jennings Report** for a different data range, click the **calendar icon** in the **Date Range** tool to select the desired beginning and end dates.



**Note:** By default, the range of the **Levey-Jennings Report** begins with the first Baseline and ends with the most recent Performance Test available for the selected Baseline.

If the default range exceeds 180 days, then the report shows data for the most recent 180-day period.

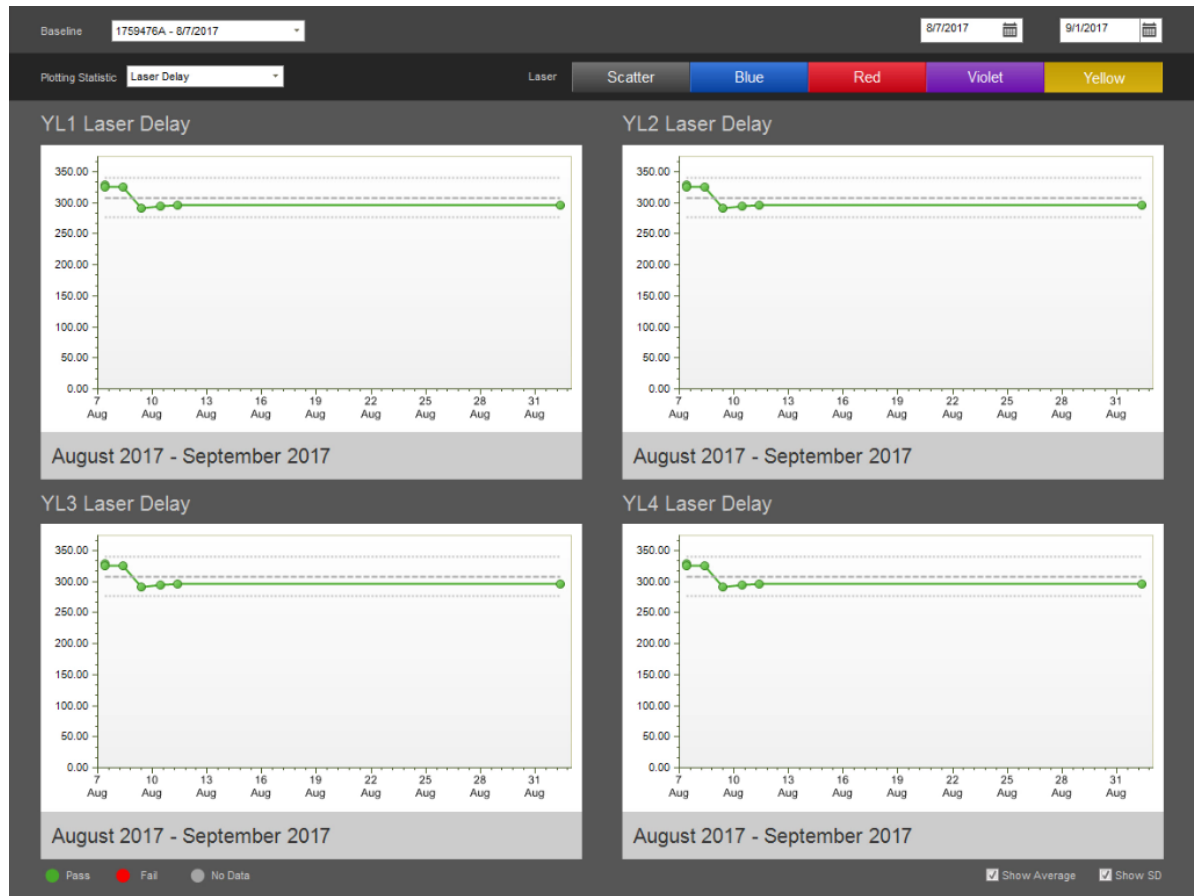
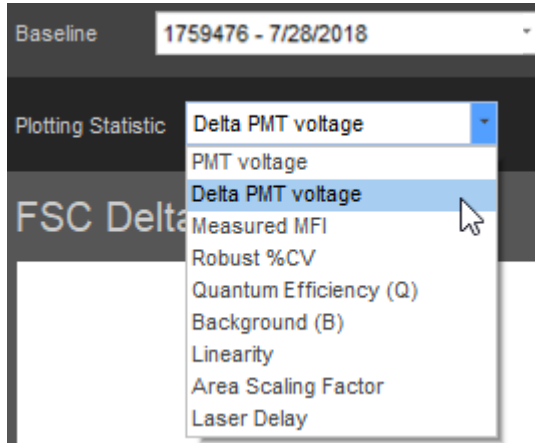
Using the **Date Range** tool, you can select a date range of up to 180 days.

If more than one result is available on a single date and the range is set such that the results overlap, the most recent result is displayed first.

## Levey-Jennings plots

- The Attune™ Cytometric Software generates **Levey-Jennings plots** for the parameters listed in Chapter 20, “Performance Test reports”.

Select the parameter of interest from the **Plotting Statistic** dropdown list to view the **Levey-Jennings plot** for that parameter.



- The **Laser** selection tool on the **Levey-Jennings Report** screen allows data to be shown for the scatter channels, blue channels, red channels, violet channels, and yellow channels.

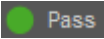

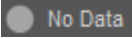
- The buttons on the **Laser** selection tool are color-coded based on the laser color. To view the **Levey-Jennings plot** for the scatter/laser channel of interest, click the button for the desired channel.



**Note:** The software shows only the channels that were configured at the time that the Performance Test was run.

- The Levey-Jennings plots** show the **Pass/Fail status** for the selected parameter within the selected date range. Each data point on the plot corresponds to a single date and the color of the data point indicates its status (see below).

If more than one result is available on a single date and the range is set such that the results overlap, the most recent result is displayed first.

Icon	Pass/Fail status
	<b>Pass:</b> Performance of the selected channel for the selected parameter is within the limits set by the Baseline calculation.
	<b>Fail:</b> Performance of the selected channel for the selected parameter deviates from the target set by the Baseline calculation.
	<b>No data:</b> There is no data for the selected channel or the parameter on that date.

- Place the pointer over a data point to view the corresponding data and the date. The decimal places are only shown for Robust %CV, Qr, Linearity, and Area Scaling Factor (2 decimal places for r%CV and Area Scaling Factor, and 3 decimal places for Q and Linearity).
- Click a data point within the **Levey-Jennings plot** to open the **Baseline** or **Performance Test result** for that date, depending on whether the data point represents the Baseline or the Performance Test.
- Use the mouse wheel to zoom in on the X-axis, which displays a scroll bar when the zoom function is applied. The Y-axis cannot be zoomed.
- If the date range is changed, the zoom is reset to the default.
- Select **Show Average** below the **Levey-Jennings plot** to view the average for the selected parameter and channel within the selected date range. The average appears as a dashed line.

Show Average

- Select **Show SD** (standard deviation) to show the standard deviation for the data. Two standard deviations appear as dotted lines above and below the center line.

Show SD

## Print the report

Select **File** ▶ **Print** to print the report.

The printout of the **Levey-Jennings Report** includes a header containing the date range of the report and the instrument information, and the Levey-Jennings plots for all configured lasers.

The first page of the printout contains both scatter channels and the first laser channel. The remaining channels are printed individually on the subsequent pages.

The following examples show the first and second pages of a **Levey-Jennings Report** printout.

Attune™ NxT Cytometer Levey-Jennings Report

3/9/2018 -7/20/2018

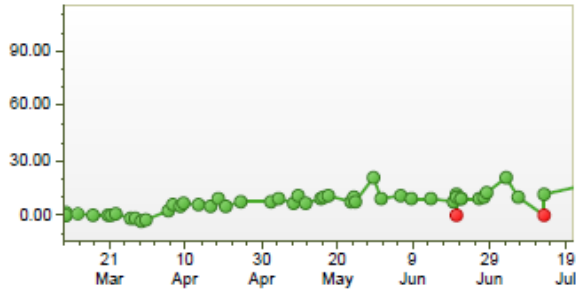
Cytometer: Attune NxT Acoustic Focusing Cytometer (Lasers: BRVY)  
Instrument Model Number: 4486521  
Instrument Serial Number: AFC2\_INVALID\_SERIAL

User: admin  
Institution:  
Software: 2.6.580.1

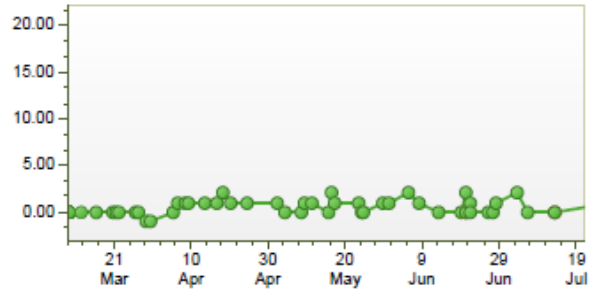
● Pass ● Fail ● No Data

Scatter Channels:

FSC Delta PMT voltage



SSC Delta PMT voltage





Attune™ NxT Cytometer Levey-Jennings Report

3/9/2018 -7/20/2018

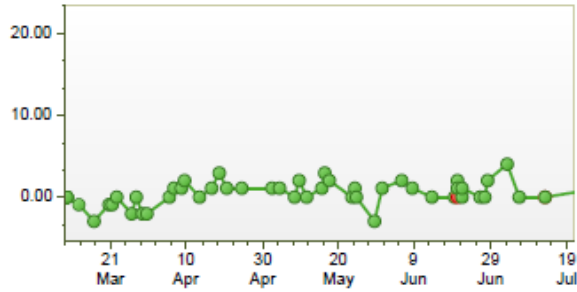
Cytometer: Attune NxT Acoustic Focusing Cytometer (Lasers: BRVY)  
Instrument Model Number: 4486521  
Instrument Serial Number: AFC2\_INVALID\_SERIAL

User: admin  
Institution:  
Software: 2.6.580.1

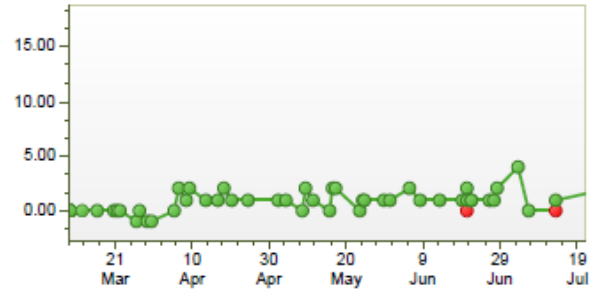
● Pass ● Fail ● No Data

Blue Laser Channels:

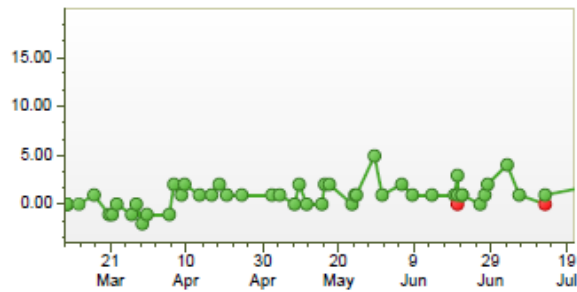
BL1 Delta PMT voltage



BL2 Delta PMT voltage



BL3 Delta PMT voltage



## Baseline Report

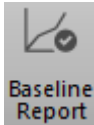
The **Baseline Report** shows the results of a selected Baseline.

When establishing a Baseline, the Attune™ Cytometric Software generates data for the parameters listed in Chapter 20, “Performance Test reports”, to determine the initial status of the cytometer and to suggest optimal minimum voltage settings for the instrument. The results of the Baseline are then used as a benchmark when tracking the daily performance of the instrument with Performance Tests.

### View the Baseline Test Results

The **Baseline Test Results** screen automatically appears after a Baseline is completed.

Alternatively, click **Baseline Report** on the **Performance Test** tab of the **BL/PT module** (“Baseline and Performance Test (BL/PT) module” on page 541).



**Baseline Test Results**

✓ Baseline test successful


Baseline: 2029773 - 1/30/2024 1/30/2024 5:26:33 PM

Channel	PMTV	Optimal PMTV	Target MFI	MFI	Robust %CV	Qr	Background	Linearity	ASF	Laser Delay	Result
FSC	369	--	300,000	298,705	2.24 %	0.000	0	0.000	1.00	1100	✓
SSC	275	--	300,000	301,083	4.42 %	0.000	0	0.000	1.00	1100	✓
BL1	420	403	300,000	295,946	1.63 %	0.037	169	0.999	1.00	1100	✓
BL2	378	394	300,000	295,983	1.50 %	0.042	173	1.000	1.00	1100	✓
BL3	457	461	300,000	300,682	2.97 %	0.015	76	1.000	1.00	1100	✓
RL1	355	367	300,000	297,420	2.98 %	0.076	44	0.996	0.97	1539	✓
RL2	381	382	300,000	300,970	2.97 %	0.011	295	0.996	0.97	1539	✓
RL3	434	410	300,000	297,070	3.22 %	0.046	124	1.000	0.97	1539	✓
VL1	249	289	300,000	300,818	0.83 %	0.019	3067	1.000	1.10	701	✓
VL2	291	293	300,000	295,296	1.01 %	0.015	1033	0.997	1.10	701	✓
VL3	373	416	300,000	297,070	1.48 %	0.018	93	1.000	1.10	701	✓
VL4	398	414	300,000	299,522	2.65 %	0.002	283	0.990	1.10	701	✓
YL1	436	438	300,000	297,844	1.52 %	0.109	59	1.000	0.91	313	✓
YL2	410	423	300,000	296,356	2.50 %	0.068	38	1.000	0.91	313	✓
YL3	468	456	300,000	296,884	3.09 %	0.013	197	1.000	0.91	313	✓
YL4	475	479	300,000	297,090	3.58 %	0.005	369	1.000	0.91	313	✓

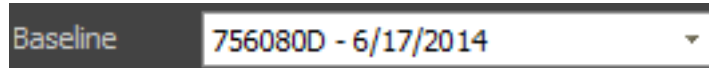
## Baseline Results

- The overall **Pass/Fail status** of the current Baseline is displayed by the **Pass/Fail indicator**, which includes a message and an indicator icon.

**Pass:**  Baseline test successful Data from all channels meet the criteria specified by the instrument.

**Fail:**  Baseline test completed with errors Data from one or more of the channels fail to meet the criteria specified by the instrument.

- To view the report for another Baseline run and Bead lot, select the Baseline of interest from the **Baseline** dropdown list.



Each **Baseline** in the dropdown list is listed in the following format:

**###** = Performance Test bead lot number

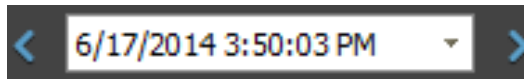
**DD/MM/YYYY** = The date Baseline status was established

---

**Note:** The **Baseline** dropdown list is active only if more than one Baseline has been established. The Baselines are ordered chronologically, with the most recent displayed first.

---

- To view results from other Baseline calculations, select the desired test date from the **Baseline** dropdown, or click the blue **arrows** to navigate through the **Baseline Reports**.



---

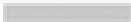
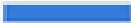




**Note:** If there are multiple reports for a single day, the reports are displayed chronologically, with the most recent displayed first.

---

## Baseline Results table

The **Baseline Results table** shows all statistics and calculations for all configured channels on the instrument.

Channel	PMTV	Optimal PMTV	Target MFI	MFI	Robust %CV	Qr	Background	Linearity	ASF	Laser Delay	Result
FSC	369	--	300,000	298,705	2.24 %	0.000	0	0.000	1.00	1100	✓
SSC	275	--	300,000	301,083	4.42 %	0.000	0	0.000	1.00	1100	✓
BL1	420	403	300,000	295,946	1.63 %	0.037	169	0.999	1.00	1100	✓
BL2	378	394	300,000	295,983	1.50 %	0.042	173	1.000	1.00	1100	✓
BL3	457	461	300,000	300,682	2.97 %	0.015	76	1.000	1.00	1100	✓
RL1	355	367	300,000	297,420	2.98 %	0.076	44	0.996	0.97	1539	✓
RL2	381	382	300,000	300,970	2.97 %	0.011	295	0.996	0.97	1539	✓
RL3	434	410	300,000	297,070	3.22 %	0.046	124	1.000	0.97	1539	✓
VL1	249	289	300,000	300,818	0.83 %	0.019	3067	1.000	1.10	701	✓
VL2	291	293	300,000	295,296	1.01 %	0.015	1033	0.997	1.10	701	✓
VL3	373	416	300,000	297,070	1.48 %	0.018	93	1.000	1.10	701	✓
VL4	398	414	300,000	299,522	2.65 %	0.002	283	0.990	1.10	701	✓
YL1	436	438	300,000	297,844	1.52 %	0.109	59	1.000	0.91	313	✓
YL2	410	423	300,000	296,356	2.50 %	0.068	38	1.000	0.91	313	✓
YL3	468	456	300,000	296,884	3.09 %	0.013	197	1.000	0.91	313	✓
YL4	475	479	300,000	297,090	3.58 %	0.005	369	1.000	0.91	313	✓

- The configured channels are listed in the following order and color-coded based on the laser color.
  -  : Scatter channels (FSC and SSC)
  -  : Blue channels
  -  : Red channels
  -  : Violet channels
  -  : Yellow channels
  -  : Green channels
- The data for the following parameters is displayed for all configured channels in the order specified below.
  - PMTV (photomultiplier tube voltage)
  - Optimal PMTV
  - Target MFI (target median fluorescence intensity)
  - Measured MFI
  - Robust %CV (robust percent coefficient of variation)
  - Quantum Efficiency (Qr)
  - Background
  - Linearity
  - Area Scaling Factor (ASF)
  - Laser Delay

- The decimal places are only shown for **Robust %CV, Qr, Linearity, and Area Scaling Factor** (2 decimal places for Robust %CV and Area Scaling Factor, and 3 decimal places for Qr and Linearity).
- **Optimal PMTV** column displays the optimal voltages only for fluorescent parameters. **The software does not calculate or change values for forward scatter (FSC) or side scatter (SSC), because those values are more shape and size dependent than stain dependent.**
- A red value with a tilde (for example, ~394 in red) in the **Optimal PMTV** column means that some voltage walking steps in BFR were successful, but other steps were inconclusive. Caution is urged. A red N/A in the **Optimal PMTV** column indicates that the optimal PMTV value could not be calculated for the detector.
- You can adopt the optimal PMTV values for all detectors, including those with tilde values, to ensure optimal performance. The BFR process does not fail even if one or more detectors yields an N/A or a tilde value.
- The **Pass/Fail status** for each configured channel is indicated by the following **status indicator icons**:
  - ✓ **Pass**: All statistics and calculations meet the criteria set by the instrument.
  - ⚠ **Fail**: One or more of the calculations or statistics deviate from the criteria set by the instrument.

## Print the Baseline Results report

Select **File ▶ Print** to print the Baseline Results report.

In addition to the Baseline Results table, the printout includes a header containing the date range of the report and the instrument information.

Attune™ NxT Cytometer Baseline Report

7/28/2018 8:28:00 PM

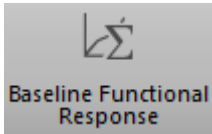
Cytometer: Attune NxT Acoustic Focusing Cytometer (Lasers: BRVY) User: admin
Instrument Model Number: 4486521 Institution: Thermo Fisher Scientific
Instrument Serial Number: 1AFC200010913 Software: 2.7.873.0

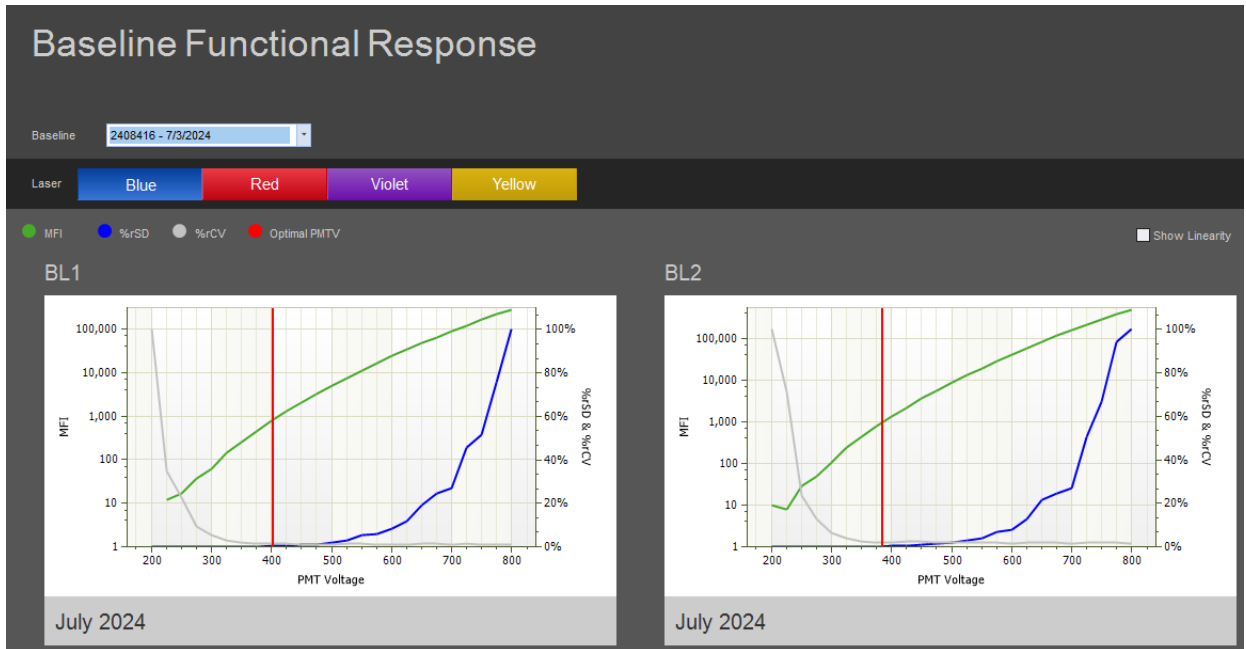
## Baseline Functional Response (BFR) Report

**Baseline Functional Response (BFR) Report** shows the BFR calculations for each fluorescent parameter in graphical form. The **BFR Report** is useful in fine tuning optimal PMT voltages and for troubleshooting inconclusive BFR results.

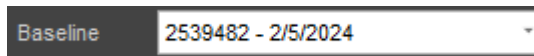
### View the Baseline Functional Response (BFR) Report

Click **Baseline Functional Response** on the **Performance Test** tab of the **BL/PT module** (“Baseline and Performance Test (BL/PT) module” on page 541).





- **Baseline** displays the date of the Baseline run and the Bead lot of the Attune™ Performance Tracking Beads used for the Baseline.  
To view the BFR report and graphs for another Baseline run and Bead lot, select the Baseline of interest from the **Baseline** dropdown list.



Each **Baseline** in the dropdown list is listed in the following format:

**###** = Performance Test bead lot number

**DD/MM/YYYY** = The date Baseline status was established

---

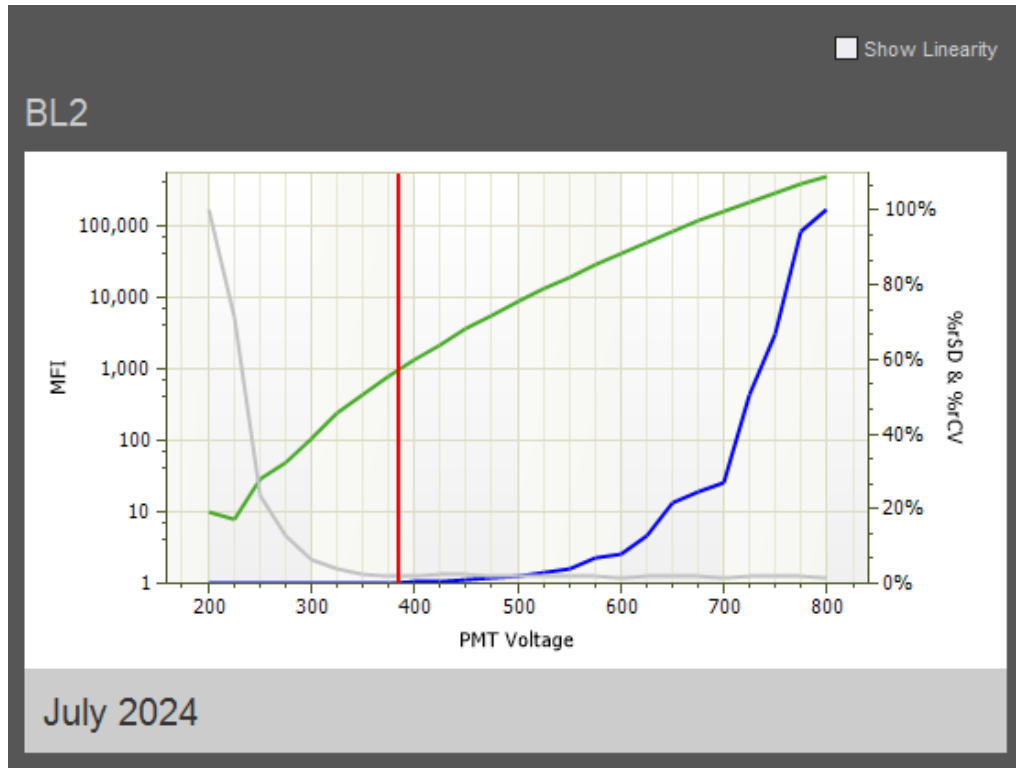
**Note:** The **Baseline** dropdown list is active only if more than one Baseline has been established. The Baselines are ordered chronologically, with the most recent displayed first.

---

- **Laser** buttons enable you select the laser for which to view the BFR results.



- **BFR plot** shows the Baseline Functional Response measurements for each fluorescent detector.



The plot displays the MFI (median fluorescence intensity), %rSD (percent robust standard deviation), r%CV (robust percent coefficient of variation) measurements and the optimum PMT voltage for the channel calculated from the BFR data (see **BFR plot legend** below).



**Note:** The BFR sets the optimal minimum PMT voltage for each channel where the robust standard deviation (rSD) of the dim bead is 3 times the rSD of the electronic noise (EN) (shown as the red line in the BFR plot).

- To show the system linearity across the PMTV range in the plot, select **Show Linearity**.

Show Linearity

**Note:** Linearity measurements provide guidance for the effective linear working range of the PMTV response. If you need to manually adjust the PMT voltages further, you can place the positive populations below the top of the linear range at the high end of the measurement.

## Print the Baseline Functional Response (BFR) report

Select **File** ► **Print** to print the BFR report.



## Overview

### Compensation

Fluorophores emit light over a range of wavelengths. Although optical filters limit the range of frequencies measured by a given detector, when two or more fluorophores are used in an experiment, there is often an overlap between the wavelength ranges.

*Compensation* is the mathematical method used to correct the emission overlap from one fluorophore into the emission channel of another fluorophore.

The Attune™ Cytometric Software Cytometric Software calculates the Compensation settings automatically as it guides you through the process. This chapter details:

- Process of setting up Compensation (“Compensation setup dialog” on page 581)
- Working with the Compensation Workspaces (“Compensation control source – wells” on page 585)
- Modifying Compensation (“Matrix dialog” on page 599)

### Compensation controls

- If the experiment requires compensation, prepare the necessary Compensation controls. You need single-stained controls (i.e., compensation beads or cells) for each fluorophore you are using for compensation. Unless you select to use a negative gate or none, you also need an unstained or isotype-labeled control.
- You can use Tubes or Plates as the source of Compensation controls. The Compensation Setup dialog provides different options for setup based on the selection of the Compensation Source. For more information, see “Compensation setup dialog” on page 581.

---

**IMPORTANT!** When you click **Record** on any of the Compensation controls in the Compensation Setup, the Instrument Settings for all fluorescent channels are shaded gray and cannot be changed. It is critical that you optimize voltages before recording any Samples or Compensation controls.

---

# Compensation setup

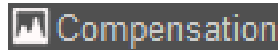
## Initiate compensation setup

To start Compensation Setup, do one of the following:

- Click the **Compensation Setup** button on the **Compensation ribbon tab** (“View tab” on page 77).



- Double-click the **Compensation node** in an Experiment on the **Experiment Explorer** (Chapter 11, “Experiment Explorer”) when there are no Compensation controls present.



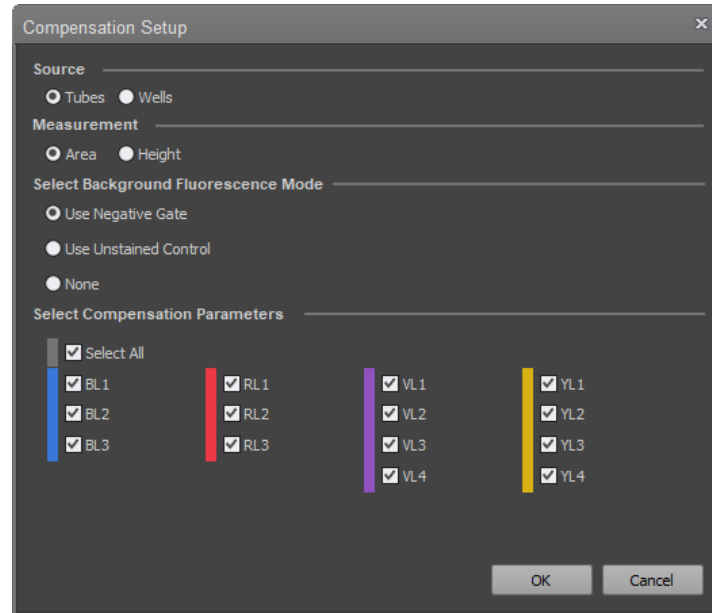
- Click the **New Compensation Control** button on the **Experiment ribbon tab** to define the Compensation samples.



Each of these methods launches the **Compensation Setup dialog**.

## Compensation setup dialog

The *Compensation Setup dialog* enables you to select the Compensation Source, the Compensation Measurement, the Background Fluorescence mode, and the required Compensation Parameters. This dialog also enables you to make modifications to an existing Compensation Setup.

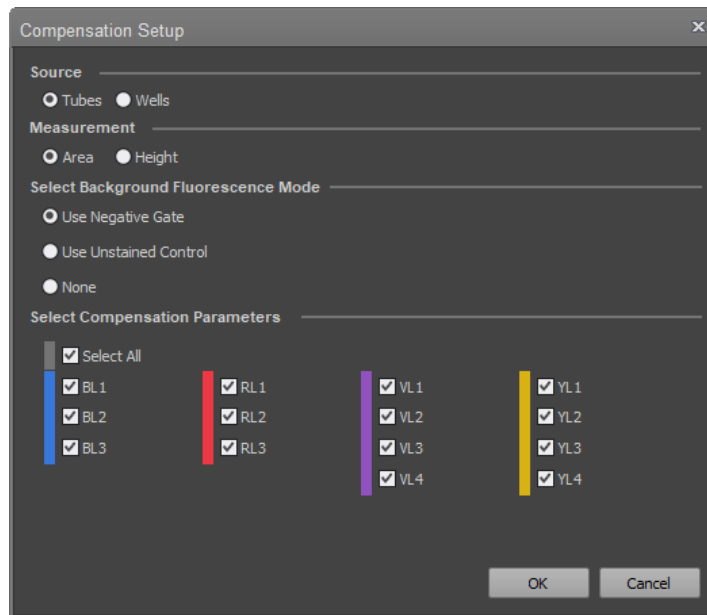


- The **Source** section shows the option for the source of Compensation controls, which is **Tubes** (“Compensation control source – Tubes” on page 581).
- The options and controls available in the remaining sections of the dialog (**Measurement**, **Background Fluorescence Mode**, and **Compensation Parameters**) are based on the Source selection.
- After Compensation controls have been created and recorded, the Source selection cannot be changed and this option becomes disabled.

## Compensation control source – Tubes

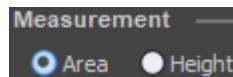
When **Tubes** are selected as the **Source** for compensation controls, the Compensation Setup dialog appears and contains the following controls:

- **Measurement** buttons (“Compensation control source – Tubes” on page 581)
- **Select Background Fluorescence Mode** buttons (“Select background fluorescence mode” on page 586)
- **Select Compensation Parameters** checkboxes (“Select compensation parameters” on page 588)



## Measurement

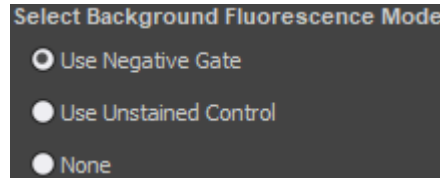
**Measurement** allows you select **Area** or **Height** as the parameter for calculating the compensation.



- By default, **Area** is selected. If no area measurements are selected in the Experiment Instrument Settings, the selection defaults to **Height**.
- When **Area** is selected, the resulting Compensation Workspace and controls use the area measurement as the parameter for calculating the compensation.  
When **Height** is selected, the resulting Compensation Workspace and controls use the height measurement as the parameter for calculating the compensation.
- If all the channels have both the height and area measurements enabled, the Measurement selection can be changed after the Compensation controls have been recorded; otherwise this option is disabled.
- If the Measurement selection is changed, the Workspaces are updated to use the new Measurement parameter (if the parameters are enabled) and the Compensation is recalculated.

## Select background fluorescence mode

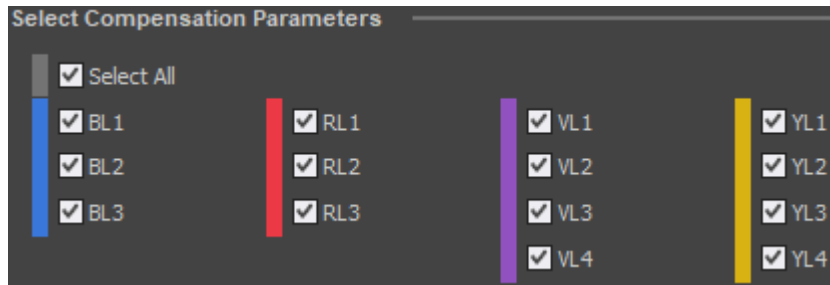
**Select Background Fluorescence Mode** enables you to select the source of the background fluorescence when calculating compensation.



- The available options are **Use Negative Gate**, **Use Unstained Control**, or **None**.
- By default, the **Use Negative Gate** is selected.
- **Use Unstained Control** adds Unstained Control as an extra Compensation control to the Compensation node in the Experiment.
- **Use Negative Gate** provides an extra Histogram gate or Rectangle gate on the Compensation Workspace for defining the negative population.
- When **None** is selected, compensation is calculated without correcting for background auto-fluorescence.
- If you change the Background Fluorescence Mode before recording any Compensation controls, the Compensation Workspaces are refreshed with the new background option and any changes are reset to the defaults.
- The Background Fluorescence Mode options are disabled after Compensation controls have been created and any Compensation controls have recorded data.  
If there is no recorded data, changing the Background Fluorescence Mode updates the Compensation controls and Workspaces accordingly.

## Select compensation parameters

**Select Compensation Parameters** contains the checkbox controls that determine the parameters used for calculating compensation.



- The **Compensation Parameter** controls are visible and enabled only for the parameters that are selected in the Parameters section of the Instrument Settings Panel, as described on “Parameters” on page 385.
- The checkboxes are arranged in columns grouped by laser color (Blue, Red, Violet, and Yellow) with a checkbox for each fluorescence detector assigned to each laser.
- By default, all enabled checkboxes are selected.
- When the **Select All** is selected, all enabled parameters are checked. When the **Select All** is deselected, all checkboxes are unchecked. Deselecting one or more of the parameters also deselects the **Select All** checkbox.
- You must select at least two Compensation Parameters. If less than two parameters are checked, the **OK** button is grayed out and a warning stating “At least two fluorescence parameters must be selected.” is shown under the buttons.



- Selection of Compensation Parameters can be modified after controls have been created and recorded. Any controls that do not have data recorded can be deselected, which removes them from the Experiment Explorer. Additional controls that have not been assigned can be selected, which adds them to the Experiment Explorer.
- The checkboxes become disabled and remain checked for any controls that have recorded data.

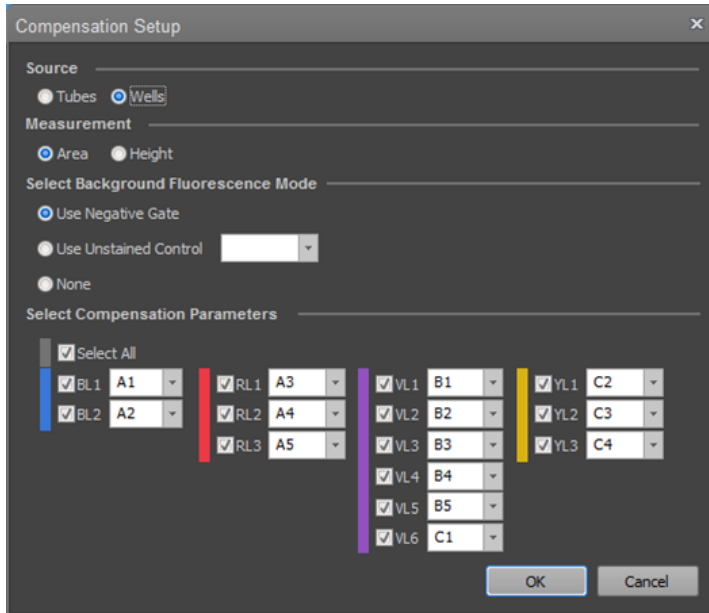
## OK and cancel buttons

- Click **OK** to close the dialog box and create or update the Compensation control for each selected parameter within the Compensation node of the Experiment in the Experiment Explorer. The Compensation Workspace for the first control or the unstained control (if **Use Unstained Control** was selected) opens automatically. An *Identity Compensation matrix*, a matrix of  $N \times N$  dimensions with ones on the main diagonal and zeros elsewhere, is created, where  $N$  is the number of compensation parameters.
- Click **Cancel** to close the dialog without making any changes.

## Compensation control source – wells

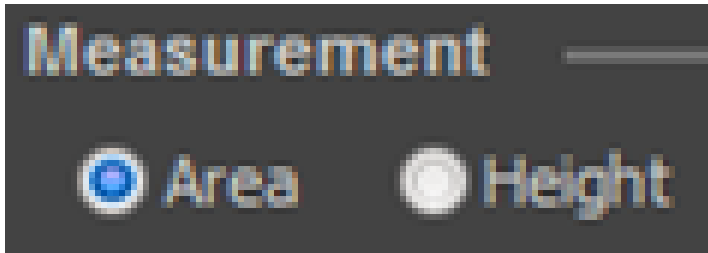
When **Wells** is selected as the Compensation control source, the Compensation Setup dialog appears and contains the following controls:

- *Measurement* radio buttons: Height, Area (“Compensation control source – wells” on page 585)
- *Select Background Fluorescence Mode* radio buttons: Negative gate, Unstained control, None (“Select background fluorescence mode” on page 583)
- *Select Compensation Parameters* combo boxes (“Select compensation parameters” on page 584)



## Measurement

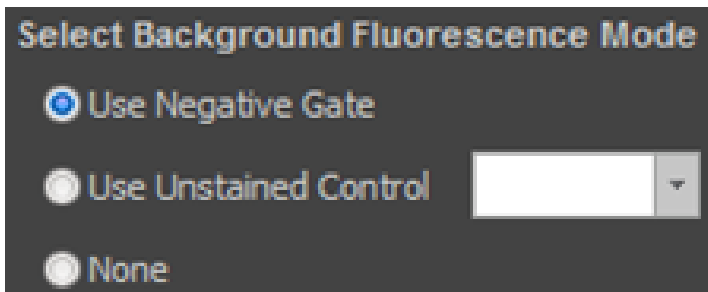
**Measurement** allows you select **Area** or **Height** as the parameter for calculating the compensation.



- By default, the **Area** is selected. If no area measurements are selected in the Experiment Instrument Settings, the selection defaults to Height.
- When **Area** is selected, the resulting Compensation Workspace and controls use the area measurement as the parameter for calculating the compensation. When **Height** is selected, the resulting Compensation Workspace and controls use the height measurement as the parameter for calculating the compensation.
- If all the channels have both the height and area measurements enabled, the Measurement selection can be changed after the Compensation controls have been recorded; otherwise this option is disabled.
- If the Measurement selection is changed, the Workspaces update to use the new Measurement parameter (if the parameters are enabled) and the Compensation is recalculated.

## Select background fluorescence mode

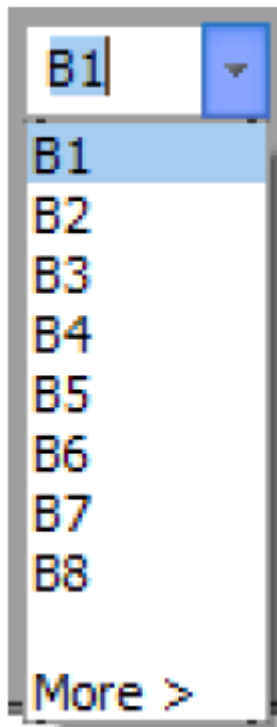
**Select Background Fluorescence Mode** allows you to choose the source of the background fluorescence when calculating compensation.



- The available options are **Use Negative Gate**, **Use Unstained Control**, or **None**. By default, the **Use Negative Gate** is selected.
- **Use Unstained Control** adds Unstained Control as an additional Compensation control to the Compensation node in the Experiment.



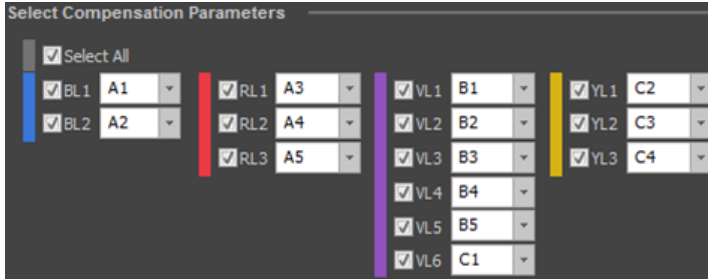
- The **Unstained Control combo box** in the Select Compensation Parameters section displays a list of all selected wells on the Heat Map and a More option to expand the list to include all wells in the plate that do not contain data.



- The wells in the combo box dropdown are listed in the same order as the Plate run order (A1–A12, B1–B12, etc.). By default, the first selected well is set as the Unstained Control
- **Use Negative Gate** provides an additional Histogram gate or Rectangle gate on the Compensation Workspace for defining the negative population (see “Compensation workspace” on page 590).
- When **None** is selected, compensation is calculated without correcting for background auto-fluorescence.
- The Background Fluorescence Mode options are disabled once Compensation controls have been created and any Compensation controls have data recorded.  
 If data has not been recorded, changing the Background Fluorescence Mode option updates the Compensation controls and compensation Workspaces accordingly.

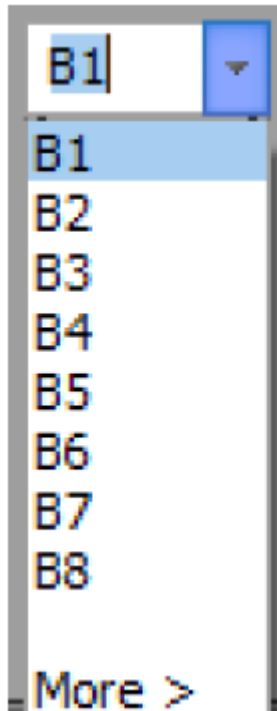
## Select compensation parameters

**Select Compensation Parameters** contains checkbox controls that determine the parameters used for calculating compensation.



- The **Compensation Parameter** controls are visible and enabled only for the parameters that are selected in the Parameters section of the Instrument Settings Panel (“Parameters” on page 385).
- The checkboxes are arranged in columns grouped by laser color (Blue, Red, Violet, and Yellow) with a checkbox for each fluorescence detector assigned to each laser.
- The checkbox labels use the display name set in the Options dialog (“Naming Options” on page 644). The display name can be modified to display the Parameter name (e.g., BL1-A), the Fluorophore name (e.g., FITC), or both (e.g., BL1-A – FITC) (“Naming Options” on page 644).
- When wells are selected prior to opening the Compensation Setup dialog, the corresponding Compensation control parameters are pre-selected on the Compensation Setup dialog and mapped to the wells in the order of well location (top to bottom, column by column).
- When wells are not preselected prior to opening the compensation setup dialog, the selected Compensation control parameters are mapped to the wells in the order of well location (top to bottom, column by column).
- To manually map a Compensation Parameter to a specific well, select the well location from the dropdown menu next to the selected parameter.

- The **Compensation Parameter** combo boxes contain a list of all selected wells and a More option to expand the list to include all wells in the plate that are unmapped.



- The wells in the combo box dropdown are listed in the same order as the Plate run order (A1–A12, B1–B12, etc.). If the list shows more than 25 well locations, a scroll bar is displayed, allowing you to view the entire list.  
 If no wells are selected, the combo box list includes all unmapped wells.
- When the **Select All** checkbox is selected, all enabled parameters are checked.



When the **Select All** checkbox is deselected, all checkboxes are unchecked. Deselecting one or more of the parameters also deselects the Select All checkbox.

- At least two Compensation Parameters must be selected for compensation to be calculated and applied to an experiment.  
 If less than two parameters are checked, the OK button is grayed out and a warning stating “At least two fluorescence parameters must be selected.” is shown under the buttons.



- Selection of Compensation Parameters can be modified after controls have been created and recorded.
  - To remove Compensation controls from the experiment, open the Compensation Setup dialog and deselect the parameter by unchecking the checkbox next to the parameter name. Only controls that do not have data can be removed.
  - To add additional controls to the compensation matrix, open the Compensation Setup dialog, select the appropriate parameter from the dialog, then choose the appropriate well from the dropdown menu next to the parameter name.

You can designate only unmapped wells as Compensation control wells. The additional Compensation controls will be listed in the Experiment Explorer and the Heat Map after clicking **OK** and closing the compensation setup dialog.
- The checkboxes become disabled and remain checked for any controls that have recorded data. If the FCS file is removed, these controls will become enabled for deselection.

## OK and cancel buttons

- Clicking **OK** closes the dialog box and creates or updates the Compensation control for each selected parameter within the Compensation Node of the Experiment in the Experiment Explorer. The Compensation Workspace for the first control or the unstained control (if **Use Unstained Control** was selected) opens automatically following compensation setup. An *Identity Compensation matrix* is created before Compensation controls are recorded. This matrix is composed of all selected compensation parameters and has values of 0%, because compensation has not yet been measured and calculated.
- Clicking **Cancel** closes the dialog and any changes are reverted.

## Compensation workspace

### Overview

*Compensation Workspace* is a special Workspace that contains all the necessary plots and gates to optimize Instrument Settings for the Compensation controls, to record Compensation controls, and to define the populations used when calculating compensation.

### Customize the compensation workspace

The Compensation Workspace has a fixed landscape view where plot layout or size cannot be adjusted. The default plot attributes are as described in the *Options dialog* (“Fonts and Styles” on page 649).

- The **Workspace Zoom** tool on the *View Ribbon tab* (“View tab” on page 77) is available for the Compensation Workspace. This tool displays all plots present in the filmstrip area of the Workspace view (“Previews” on page 132) and allows you to zoom in and out of each plot.

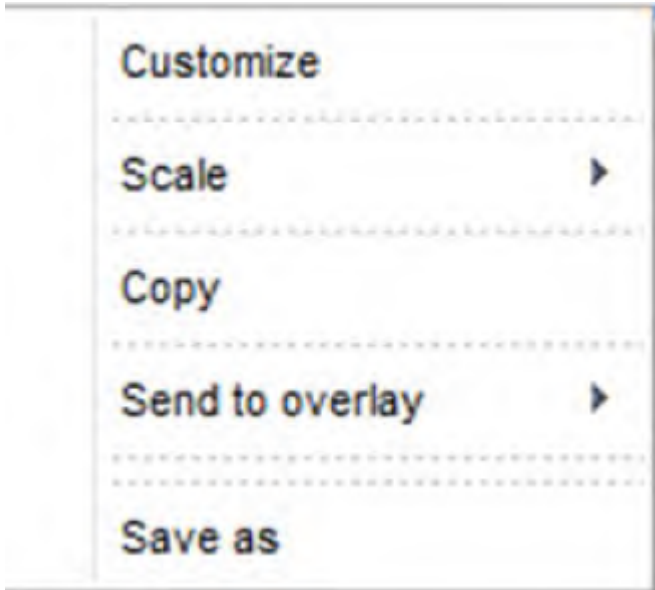




- Plot scale ranges for all parameters are set to **manual**, which allows you to enter the minimum and maximum values manually. The allowable range depends on the type of scale selected. The default maximum value for each parameter is determined by the range set in the FCS file. By default, the maximum is set to  $2^{26}$  for Event and Time counts, to  $2^{20}$  for Height and Area measurements, and to  $2^{10}$  for Width measurements. The minimum value for logarithmic parameters is set to 1, and the minimum for linear parameters is set to 0.
- The plot attributes listed below can be customized as described in “Customize Plot Options” (“Customize plot options” on page 421) and in “Workspace view” (Chapter 5, “Workspace view”).
  - Resolution
  - X and Y parameters on gating plots and where specified below
  - Scale type and range
- The plot and axis titles cannot be modified.
- The gates used on the Compensation Workspace plots use the gate style and gate label style default settings as described in the Options dialog (“Display Parameter Name as” on page 645).
- Gates cannot be customized and their names cannot be modified.

## Plot context menu

*Plot context menu* is displayed when you right-click on any empty area (i.e., white space) within the plot boundary, but not on another active area (gate, plot title, or plot axis).

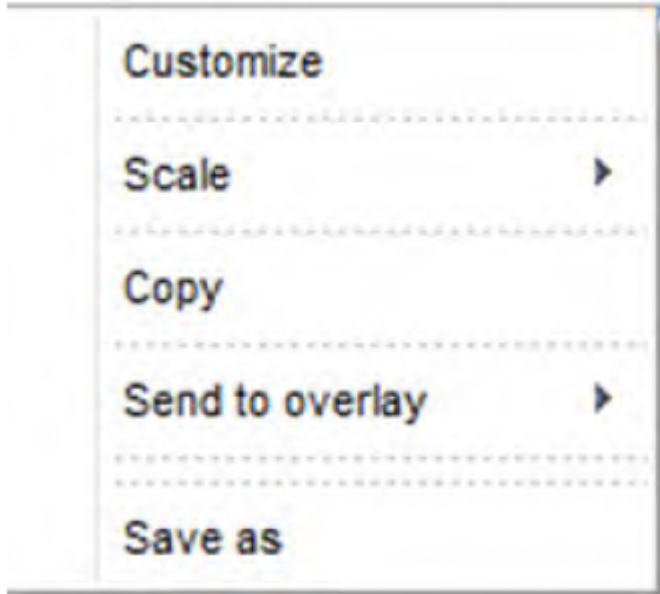


- Plot context menu contains the following options:
  - Customize
  - Scale
  - Copy
  - Send to Overlay
  - Save As
- All options listed function as described in “Plot context menu” on page 189.

## Plot axis context menu

*Plot axis context menu* is displayed when you right-click on a plot scale or parameter name. The plot axis context menu options displayed depend on the plot axis selected.

- The X-axis parameter of fluorescence plots cannot be changed, and only the following options are displayed:
  - Customize
  - Scale
  - Copy
  - Send to Overlay
  - Save As



- For the Scatter plots and Y-axis of dual-parameter fluorescence plots, the parameter can be changed, and the following options are displayed:
  - Customize
  - Scale
  - Parameter
  - Copy
  - Send to Overlay
  - Save As



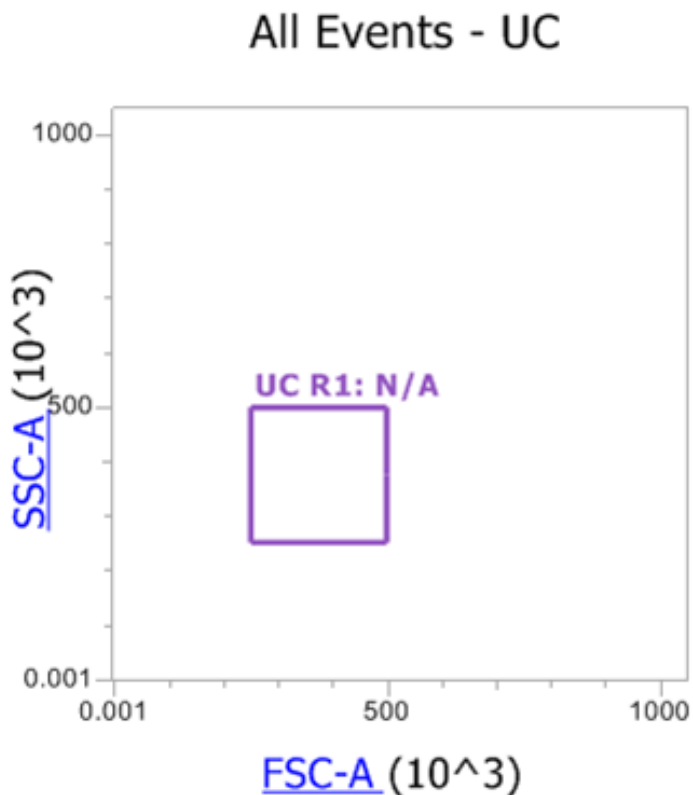
- All options listed function as described in “Plot context menu” on page 189.

## Gate context menu

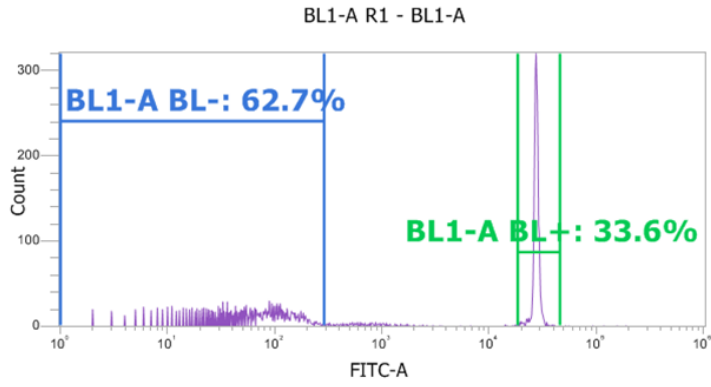
- The Gate context menu in the Compensation Workspace is available only for gates on dual-parameter scatter plots.
- Right-clicking on any part of a gate, on the gate boundary line, or a control point displays the Gate context menu.
- The context menu for dual-parameter scatter plot gates consists of a single option: Apply gate shape to all controls.
- This option applies the current scatter gate type, size, and location to the scatter gate on all Compensation controls within the current Experiment.
- Apart from the scatter gate described above, there is no gate context menu for compensation plots.

## Compensation workspace for histogram view

The Workspace for the *Histogram View* consists of a *Gating plot* and a *Histogram Compensation plot* corresponding to the selected Compensation Parameter.

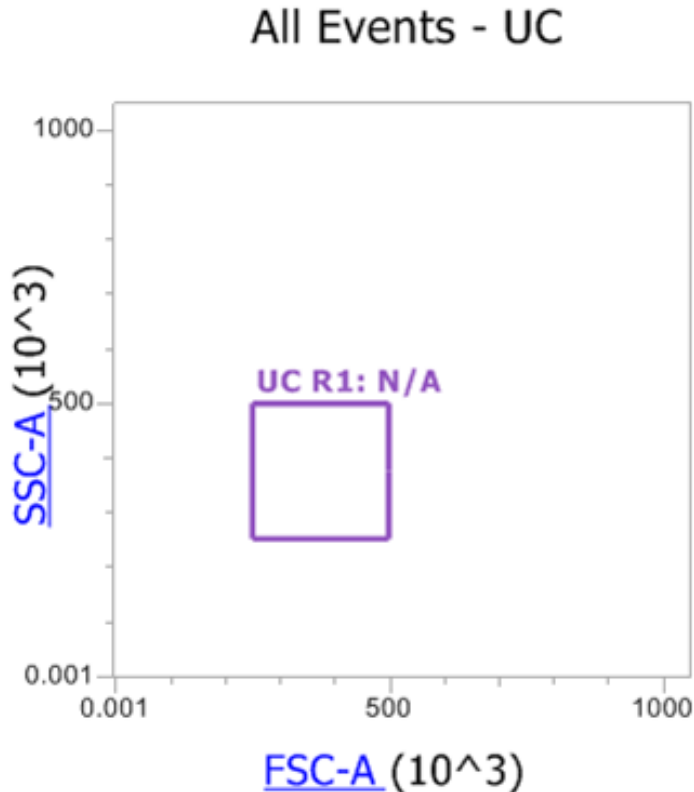






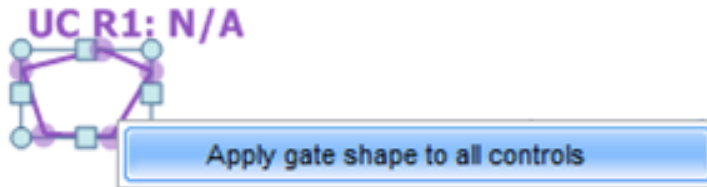
## Gating plot

The *Gating plot* is an ungated dual-parameter density plot used to gate the events shown on the Compensation plot.



- By default, the parameters for the X- and Y-axes are set to FSC and SSC, respectively, and use the measurement specified in the Measurement section of the Compensation Setup dialog (“Compensation setup dialog” on page 581).
- The Gating plot type can be changed to Ellipse or Contour Autogate.
- The Gating plot can only be set to display *All Events*. The Plot title hyperlink, if shown, is disabled.
- The Parameter names on the Gating plot are hyperlink controls that link to the Parameter selection dropdown list (see “Plot parameter hyperlink” on page 205). You can change the selected parameters for both the X- and Y-axes.
- The Gating plot contains a single 5-point polygon gate (drawn in a rectangle) named *Parent*. This gate cannot be deleted.
- The default position of the Parent gate has the following (X, Y) coordinates:
  - lower left corner coordinates (250000, 250000)
  - upper right corner coordinates (500000, 500000)
- You can move and resize the Parent gate as needed.
- Any adjustment made to the Parent gate automatically updates the Histogram plots and the calculated Compensation matrix (“Matrix dialog” on page 599).

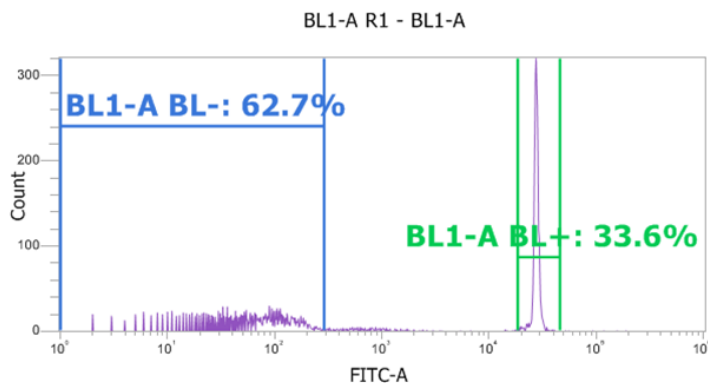
- To apply the position of the Parent gate to other compensation controls, right-click on the gate and select **Apply gate shape to all controls**.



- Initially, when the Gating plot gate position is moved on the Workspace of the first Compensation control, it automatically applies to all Compensation control samples. After a gate is moved on subsequent Compensation control samples, moving the gate on the Workspace of the first control has no effect.

## Histogram compensation plot without negative gate

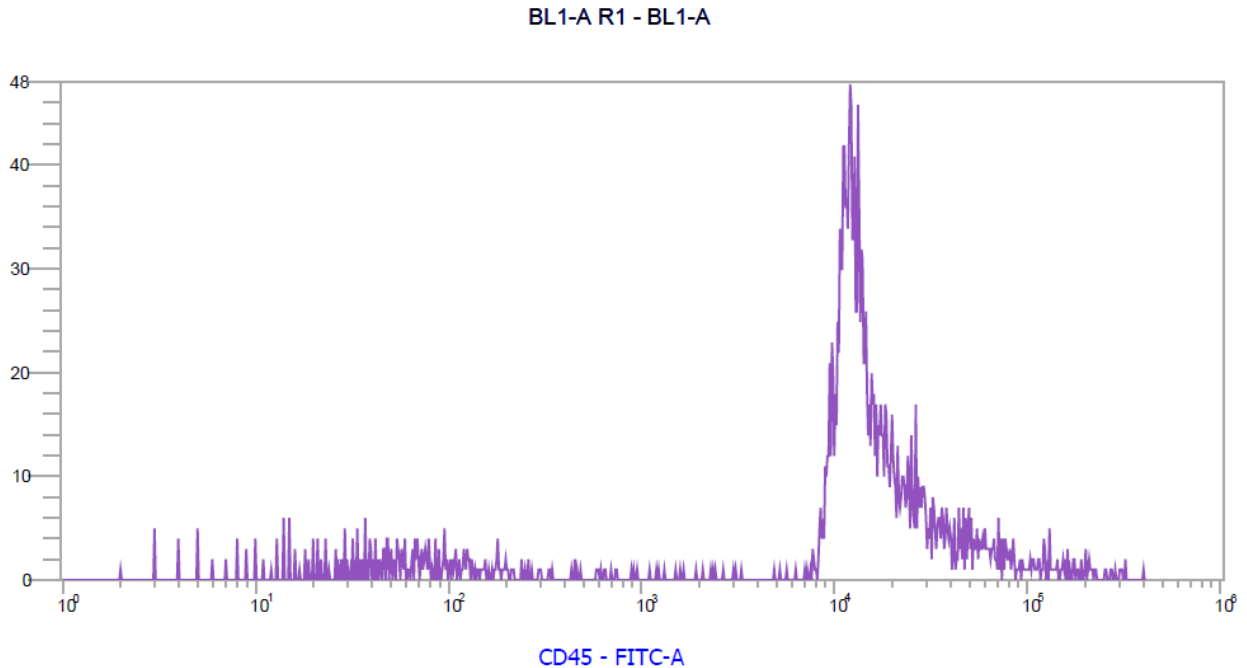
The *Histogram Compensation plot* is a single-parameter daughter plot derived from the Parent gate on the dual-parameter Gating plot (“Gating plot” on page 596). This plot type cannot be changed.



- The X-axis parameter on the plot is the currently selected Compensation control on the Experiment Explorer. This parameter cannot be changed.
- The Gate name hyperlink on the Plot title, if shown, is disabled.
- The Compensation plot contains a single Histogram gate, which is named in the following format: “<parameter channel name> +”. For example, BL1+.
- The default position of the Histogram gate has the following (X, Y) coordinates:
  - left coordinate (500,000)
  - right coordinate (750,000)
- You can move and resize the gate as needed.
- The Compensation Histogram gate cannot be deleted.
- Any adjustment made to the Histogram gate automatically updates the calculated Compensation matrix (“Matrix dialog” on page 599).

## Histogram compensation plot with negative gate

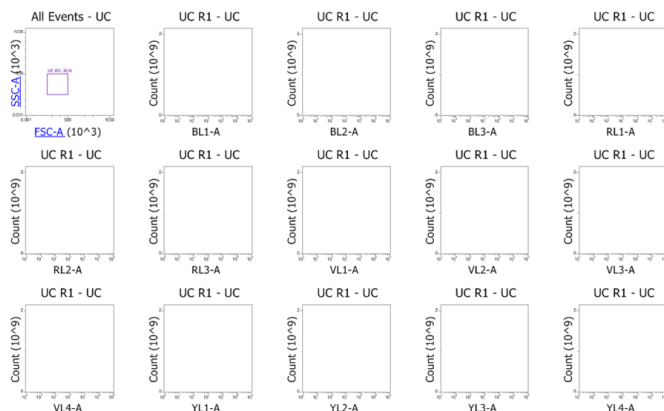
This Compensation plot is identical to the Histogram Compensation plot (“Histogram compensation plot without negative gate” on page 597) with the exception that it contains an additional gate defining a negative population.



- When **Use Negative Gate** is selected in the Compensation Setup dialog (“Compensation setup dialog” on page 581), the Histogram plot contains an additional gate to define a negative population.
- The negative gate is named in the following format: “<parameter channel name> -”. For example, BL1-.
- The default position of the Negative gate has the following (X, Y) coordinates:
  - left coordinate (1)
  - right coordinate (1000)

## Unstained control workspace

When **Use Unstained Control** is selected in the Compensation Setup dialog (“Compensation setup dialog” on page 581), an additional Control Workspace for the Unstained control is created.



- The Unstained Control Workspace contains a Gating plot as described on “Gating plot” on page 596 and Histogram plots for each Compensation parameter.
- The Histogram plots for Compensation parameters on the Unstained Control Workspace do not display any gates. All Histogram plots are gated on the Parent gate.
- The background signal is calculated based on the median of the gated data for each parameter; no additional gates are required.

## Matrix dialog

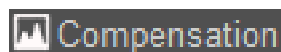
### Overview

The *Matrix dialog* shows the active compensation associated with the current Experiment or FCS file and enables you to manually edit or reset the *Spillover matrix* values.

- Click the **View Matrix** button on the *Compensation ribbon tab* to open the View Matrix dialog.



Alternatively, double-click the **Compensation node** on the Experiment Explorer when Compensation controls are present.




---

**Note:** The *Spillover matrix* shows the amount of spillover from each fluorophore into each of the other fluorescent channels.

---

View Matrix

Matrix - Experiment(9)

Spillover values are read across rows

Spillover

	BL1-A	BL2-A	RL1-A	RL2-A	RL3-A	VL1-A	VL2-A	VL3-A	VL4-A	VL5-A	VL6-A	YL1-A	YL2-A	YL3-A
BL1-A	100.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00
BL2-A	0.00	100.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00
RL1-A	0.00	0.00	100.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00
RL2-A	0.00	0.00	0.00	100.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00
RL3-A	0.00	0.00	0.00	0.00	100.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00
VL1-A	0.00	0.00	0.00	0.00	0.00	100.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00
VL2-A	0.00	0.00	0.00	0.00	0.00	0.00	100.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00
VL3-A	0.00	0.00	0.00	0.00	0.00	0.00	0.00	100.00	0.00	0.00	0.00	0.00	0.00	0.00
VL4-A	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	100.00	0.00	0.00	0.00	0.00	0.00
VL5-A	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	100.00	0.00	0.00	0.00	0.00
VL6-A	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	100.00	0.00	0.00	0.00
YL1-A	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	100.00	0.00	0.00
YL2-A	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	100.00	0.00
YL3-A	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	0.00	100.00

OK Cancel Reset

- The Matrix title indicates the source of the active compensation, which is the Experiment or the Sample name based on the selection made in the *Apply group* of the Compensation ribbon tab (“Apply group” on page 100) for the current Sample.
- If the Matrix is opened by double-clicking the **Compensation node**, the compensation source is always the Experiment.

## Matrix dialog table

The *Matrix dialog table* shows the Spillover values for the current compensation arranged in a table. The Spillover values in the table can be edited.

- The row and column labels use the Channel name (\$PnN) by default.
- The row and column labels have a background color based on the laser color associated with the channel. If a channel color is not available for a parameter (i.e., non-Attune™ file), the default fill is white and the font color is black.
- When a textbox is selected, the column and row is highlighted at the point of their intersection in blue. The diagonal values of the matrix are shaded in light gray.

	BL1-A	BL2-A	BL3-A
BL1-A	100.00	0.00	0.00
BL2-A	0.00	100.00	0.00
BL3-A	0.00	0.00	100.00
RL1-A	0.00	0.00	

- In the Spillover matrix, the textbox controls accept numbers from 0% to 100%, and allow the use of two decimal places.

The diagonal values cannot be adjusted in the Spillover matrix is shown.

- You can enter up to 7 total characters into the textboxes in the Spillover matrix, but all entered values are rounded to 2 decimal places when you click elsewhere or when you click **Enter**. Only numbers, decimals, and + and – signs can be entered in the textboxes.

- The Spillover values in the table are written across rows. For example, the spillover of FL1 into FL2 is shown in row 1, column 2.

Spillover	FL1	FL2	FL3	FL4	FL5	FL6
FL1						
FL2						
FL3						
FL4						
FL5						
FL6						

In the example below, the spillover from FL1 (BL1-A) into FL2 (BL2-A) is 38.10%.

	BL1-A	BL2-A	BL3-A	RL1-A
BL1-A	100.00	38.10	5.86	0.01
BL2-A	0.51	100.00	24.18	0.00
BL3-A	0.22	25.40	100.00	13.92
RL1-A	0.06	0.04	0.01	100.00

- When the cursor is hovered over the table textboxes, tooltips are displayed. When the Spillover matrix is displayed, the tooltip reads: “%<Row Label> in <Column Label>”. For example, %BL1 in BL2.

## OK, reset, and cancel buttons

### OK button

- OK** accepts any changes made to the matrix and applies the updated compensation values to the dataset.
- If changes are made to the matrix and **OK** is not clicked, the changes will be lost.
- If the matrix values are adjusted during recording a Sample and a stop gate is set on a population, the stop gate condition recounts the events after **OK** is clicked.

### Reset

- Reset** resets the compensation values to their default states as described below.
- If Experiment-level compensation is modified and Compensation controls are available, clicking **Reset** recalculates the compensation values based on the current Histogram gate positions on the Compensation Workspace.
- If any existing Experiment-level compensation settings are modified and Compensation controls are not available (for example, if the Experiment has been duplicated), clicking **Reset** reverts the Compensation matrix to the original Experiment-level calculated values.
- If FCS file compensation is modified, clicking **Reset** restores the compensation values to the values specified in the FCS data file.

## Cancel

- **Cancel** undoes any changes made to the Spillover matrix and retains the existing values.

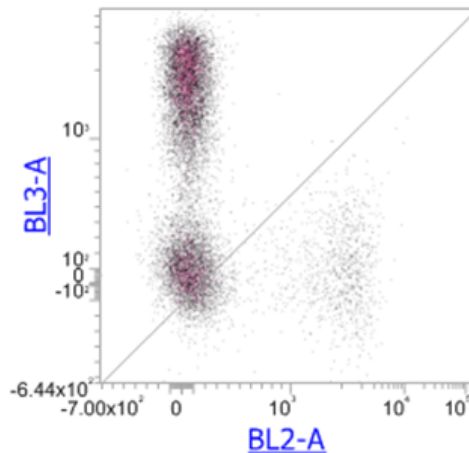
# On-Plot compensation adjustment

## Overview

- When **Plot Compensation** state button on the Compensation ribbon tab is turned **ON**, you can drag event populations displayed on dual-parameter fluorescence plots to new positions, which changes the associated Compensation matrix values.



- When **Plot Compensation** is **ON**, hovering over a dual-parameter fluorescence plot displays a tooltip showing the compensation coefficients. The coefficients displayed are the spillover values.
- For the selected plot, a diagonal line from the lower left to the upper right distinguishes the two coefficient sections.



Dragging a selected population above the line in a horizontal direction adjusts one coefficient, dragging below the line in a vertical direction adjusts the other coefficient.

- For Spillover values, the compensation coefficients can be dragged from 0% to 100%.



## Manual on-plot compensation adjustment

- Click the selected population, then drag the population to the desired location on the plot.
- The sensitivity of the dragging action depends on where the plot is clicked.
  - Clicking and dragging the population closer to the diagonal line results in larger adjustments to the spillover values.
  - Clicking and dragging farther from the diagonal line makes finer adjustments to the spillover values.
  - The exact values of the changes depend on the scale of plot and the range in which a population is moved.
- Gate labels are not visible during the drag operation.
- Only the current plot is updated during dragging. When the dragging is complete, all plots and preview plots affected by the change are updated. The compensation adjustment is only registered into the undo stack when the dragging is complete.
- If compensation values are adjusted during recording a Sample and a stop gate is set on a population, the stop gate condition recounts the events after the mouse button has been released
- You can also adjust the compensation for the selected plot using the arrow keys. The **up** and **down arrow** keys adjust the Y-axis coefficient, and the **left** and **right arrow** keys adjust the X-axis coefficient. Each arrow key changes the compensation value by 1 unit. Holding down the arrow keys enables continual adjustment while the key is pressed.
- Holding **Shift+arrow** keys change the compensation value by 0.1 units.

## Compensation acquisition workflow

### Overview

After the Tube or Well Compensation controls have been defined, they can be acquired. The expected workflow is to go through a round of Instrument Settings optimization to correctly set the voltages, thresholds, and gates, followed by a round of recording. At this point, the recorded Compensation controls are factored into the compensation calculation.

---

**IMPORTANT!** After you have recorded all Compensation controls and calculated and applied the Compensation matrix, you cannot adjust the PMT voltages for experimental data.

---

### Compensation acquisition from tubes

- Compensation setup and acquisition from Tubes is performed manually.
- You can optimize the Instrument Settings for Compensation controls using the Unstained Control Workspace or the Workspace associated with the Compensation control sample.
- The optimization process for Instrument Settings is repeated for all Compensation controls.
- After the optimization of Instrument Settings is completed, load the first Compensation control, click on its corresponding Sample tube, then record using the Collection Panel controls (“Collection controls” on page 345). Repeat the process for each Compensation control.

- While recording, gates can be moved.
  - Adjust the Polygon parent gates so that the Parent gate includes the target population (“Gating plot” on page 596).
  - Adjust the Histogram gates to include the positive and negative populations (“Histogram compensation plot without negative gate” on page 597–“Histogram compensation plot with negative gate” on page 598).
- The Compensation matrix is calculated as new Compensation controls are recorded. The calculation will only include the controls that have been recorded.

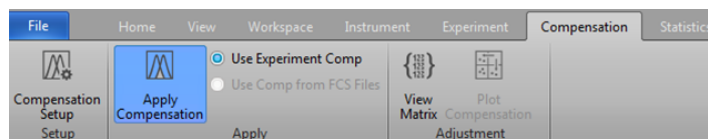
## Compensation acquisition from wells

- Compensation setup and acquisition from Wells is initiated using the **Set Up Comp** button in the **Collection Panel** (“Buttons displayed” on page 352).  
Alternatively, you can acquire Compensation controls by selecting **Record All** on the **Collection Panel** menu.
- You can navigate between Compensation controls using the **Navigation buttons** in the **Collection Panel** (“Navigation buttons” on page 355).
- You can optimize each Compensation control using the Workspace associated with the Compensation control sample.
- When **Set Up Comp** is clicked, the Sample is acquired in the Set Up mode, allowing you to adjust the voltages and thresholds.
- During the **Set Up Comp** or **Record Comp** phases, you can adjust and set any gates on the gating plot and the corresponding histogram or dual parameter plot.
- The Compensation matrix is calculated as new Compensation controls are recorded. This calculation only includes the controls that have been recorded.

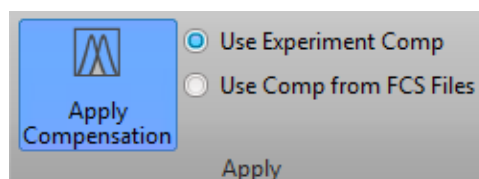
## Compensation of instrument data

For Samples acquired from the instrument, the Compensation matrix is applied to both the Area and Height measurements (if available) regardless of whether the compensation was calculated using Area or Height measurements.

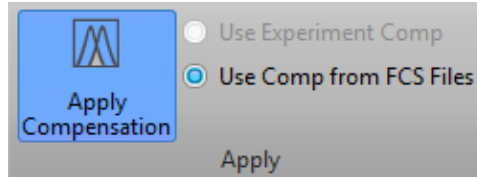
- After Compensation controls have been recorded, compensation is automatically applied to all Samples in the Experiment as indicated by the blue colored Apply Compensation button in the Compensation tab of the Ribbon bar.



- When an Experiment includes compensation and FCS files have been recorded, the **Use Experiment Comp** button becomes active.



- If FCS files have been imported into an Experiment, the compensation applied to the Sample is the compensation included in the FCS file itself. In this instance, the **Use Comp from FCS Files** button becomes active.



## New Experiment dialog

The **New Experiment dialog** enables you to create New Experiments of the following types:

- Tube-only Experiment (“Create a Tube-only experiment” on page 607)
- Plate Experiment (“Create a plate experiment” on page 609)

## Open the New Experiment dialog

On the **Ribbon bar**, click **File** to open the **File menu**, then select **New Experiment**.

Alternatively, click the **New Experiment** button in the **Home tab** or use the keyboard shortcut **Ctrl+N**.



---

**Note:** The **New Experiment** option is enabled only when the instrument is not acquiring.

---

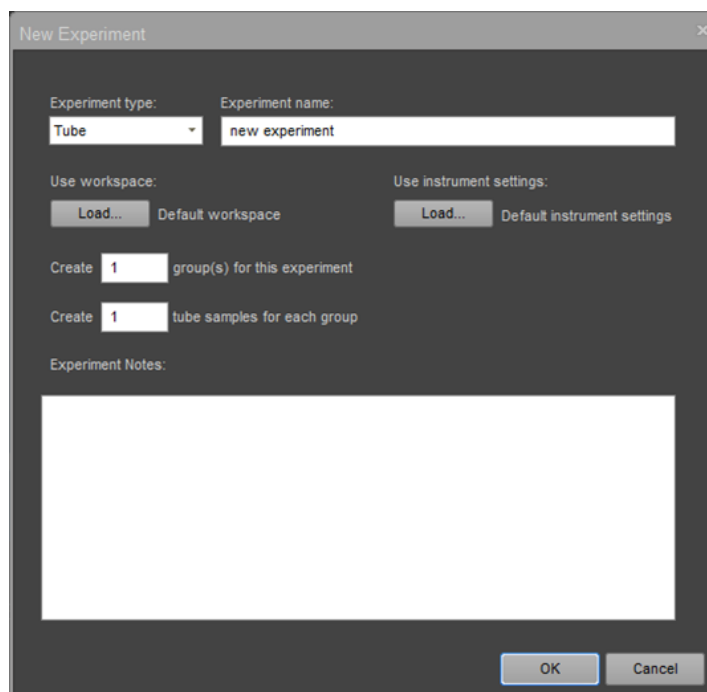


Figure 125 New Experiment dialog for Tube Experiments

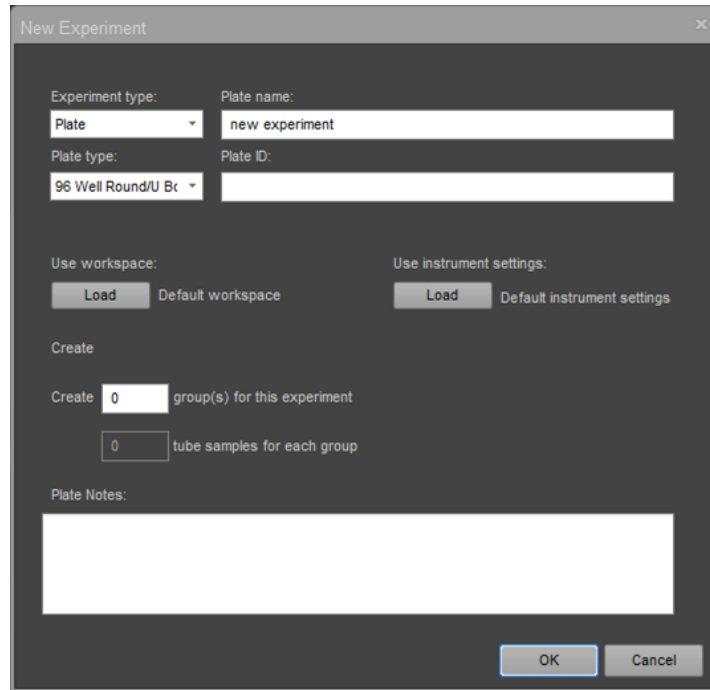


Figure 126 New Experiment dialog for Plate Experiments

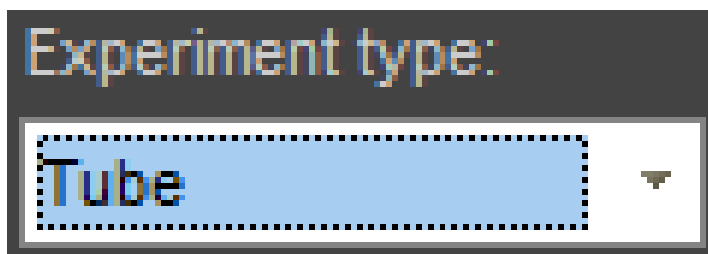
---

**Note:** The contents of the **New Experiment dialog** varies depending on the experiment type you select.

---

## Create a Tube-only experiment

1. On the **Ribbon** bar, click **File** to open the File menu, then select **New Experiment**.
2. From the Experiment type dropdown menu, select **Tube** (default selection).



The contents of the New Experiment dialog changes to display the Tube Experiment options.

3. Accept the default Experiment name or enter a new name.

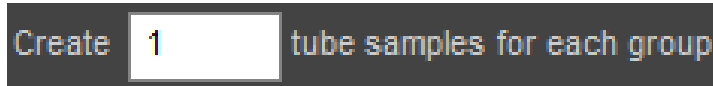
- You can enter up to 50 characters; the following characters are not permitted: \ / : \* ? " < > | % . The software automatically removes trailing or leading spaces.
- Names cannot end with a period character or be any of the following system reserved names: CON, PRN, AUX, CLOCK\$, NUL, COM1, COM2, COM3, COM4, COM5, COM6, COM7, COM8, COM9, LPT1, LPT2, LPT3, LPT4, LPT5, LPT6, LPT7, LPT8, LPT9.
- Duplicate names are not permitted.
- The default name for a Tube experiment is **Experiment**. If an Experiment using the default name already exists, a numerical suffix is added in parentheses; for example, **Experiment (2)**.

4. Accept the default **Workspace**.

5. Accept the default **Instrument Settings**.

6. Enter the number of Groups to create for the Experiment.

7. You can enter a number from **1** to **25**. The software adds the number of Groups to the Experiment Explorer.
8. The software creates a unique name for each Group with the default Group name set in the Options dialog ► General tab and appends it with a numerical suffix; for example, **Group (2)**.
9. Enter the number of Tube samples to create for each Tube group.



10. You can enter a number from **1** to **400**. If the number of Groups times the number of Samples is > 400, the number of Samples within a Group reverts to a number that is ≤400.
11. The software adds the number of Samples to Experiment Explorer.
12. The software creates a unique name for each Sample with the default Sample name set in the Options dialog ► General tab and appends it with a numerical suffix; for example, **Sample (2)**.
13. *Optional:* Enter notes for the Experiment. You can enter up to 500 characters; any character is permitted.

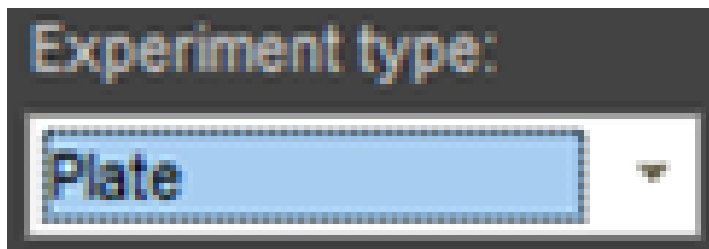


14. Click **OK** to create the New Experiment and close the New Experiment dialog. The software opens the first Sample in the Experiment and the Experiment Workspace.  
To close the dialog without creating an Experiment, click **Cancel**.

## Create a plate experiment

A *Plate Experiment* includes one Plate, and but it can also include Tube groups and Tube samples.

1. On the **Ribbon** bar, click **File** to open the File menu, then select **New Experiment**.
2. From the Experiment type dropdown menu, select **Plate**.



The contents of the New Experiment dialog changes to display the Plate Experiment options.

3. Accept the default plate name or enter a new name.

- You can enter up to 50 characters; the following characters are not permitted: \ / : \* ? " < > | % . The software automatically removes trailing or leading spaces.
  - Names cannot end with a period character or be any of the following system reserved names: CON, PRN, AUX, CLOCK\$, NUL, COM1, COM2, COM3, COM4, COM5, COM6, COM7, COM8, COM9, LPT1, LPT2, LPT3, LPT4, LPT5, LPT6, LPT7, LPT8, LPT9.
  - Duplicate names are not permitted.
  - The default name for a Plate Experiment is **Plate**. If an Experiment using the default name already exists, a numerical suffix is added in parentheses to ensure a unique name; for example, **Plate (2)**.
4. From the **Plate type dropdown** menu, select the type of plate that you are using for the Experiment.



5. Enter the plate ID.

You can enter up to 50 characters; all characters are permitted. The software automatically removes trailing or leading spaces.

6. *Optional:* If you are using a foil cover to protect the sample plate from condensation or evaporation, select **Use Foil Cover**.



- When the **Use Foil Cover** option is enabled, the autosampler disables the probe collision sensor, which allows the use of a foil cover on the plate.  
When this option is enabled, the software displays the “Foil cover present” warning dialog at the beginning of the plate acquisition to ensure that the correct plate is used.
- Use Foil Cover selection persists as part of a Template, and when exporting or duplicating a Plate.

7. *Optional:* If you are using a cooling block, select **Use Cooling Block**.



- When **Use Cooling Block** option is selected, the autosampler accounts for the extra height that the cooling block adds to the plate specification.
- When available, this option is enabled only for standard 96-well round bottom plates (such as Thermo Scientific™ 96-well Microtiter™ Microplates, Cat. No. 2205). A standard plate is defined as a plate that is within the height range of 14.35 mm ± 0.76 mm.
- Use Cooling Block selection persists as part of a Template, and when exporting or duplicating a Plate.

---

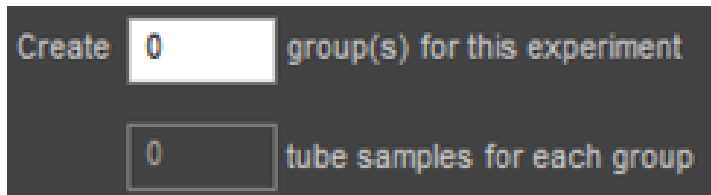
**Note:** Use Foil Cover and Use Cooling Block options are available only when a CytKick™ Max™ Autosampler is connected to the Attune™ instrument. Otherwise, they are not visible.

---

8. Accept the default **Workspace** or click **Load** to select a saved Workspace file.
9. Accept the default **Instrument Settings** or click **Load** to select a saved Instrument Settings file.



10. If you are including tubes, enter the number of **Tube Groups** to create for the Experiment.

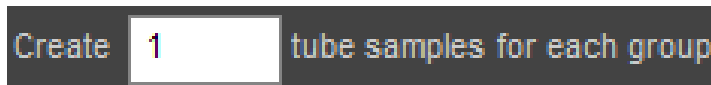


Create  group(s) for this experiment

tube samples for each group

- You can enter a number from **0** to **5**, depending on the plate type you have selected in step 4 on page 610. The default is **0**.
- The software adds the number of Groups to Experiment Explorer.
- If zero Groups are entered, then the Tube Samples field is disabled.
- The software creates a unique name for each Group with the default Group name set in the Options dialog ► General tab and appends it with a numerical suffix; for example, **Group (2)**.

11. Enter the number of **Tube Samples** to create for each Tube Group.



Create  tube samples for each group

- For a 96-well plate, you can enter a number from **1** to **304**. If the number of Groups times the number of Samples is >304, the number of Samples within a group will revert to a number that is ≤304.
- For a 384-well plate, you can enter a number from **1** to **16**. If the number of Groups time the number of Samples is >16, the number of Samples within a Group will revert to a number that is ≤16.
- The software adds the number of Samples to Experiment Explorer (“Files view” on page 287).
- The software creates a unique name for each Sample with the default Sample name set in the Options dialog ► General tab and appends it with a numerical suffix; for example, **Sample (2)**.

12. *Optional:* Enter notes for the Experiment. You can enter up to 500 characters; any character is permitted.



Notes:

13. Click **OK** to create the New Experiment and close the New Experiment dialog. The software opens the Heat Map view (Chapter 6, “Heat Map View”).

To close the dialog without creating an Experiment, click **Cancel**.

## New experiment from template dialog

The *New Experiment from Template* dialog is a modal dialog that allows you to create a new Experiment from an existing template with pre-defined settings by guiding you to select an Experiment template and enter experiment information (such Experiment name, description, notes).

The template applies the template's Workspace, Instrument settings, Run Protocols, Heat Map settings, and defined compensation to the new experiment and creates a new experiment with the same number of mapped groups and samples.

---

**Note:** For Experiments created from a template, no FCS files are attached to any of the Samples.

---

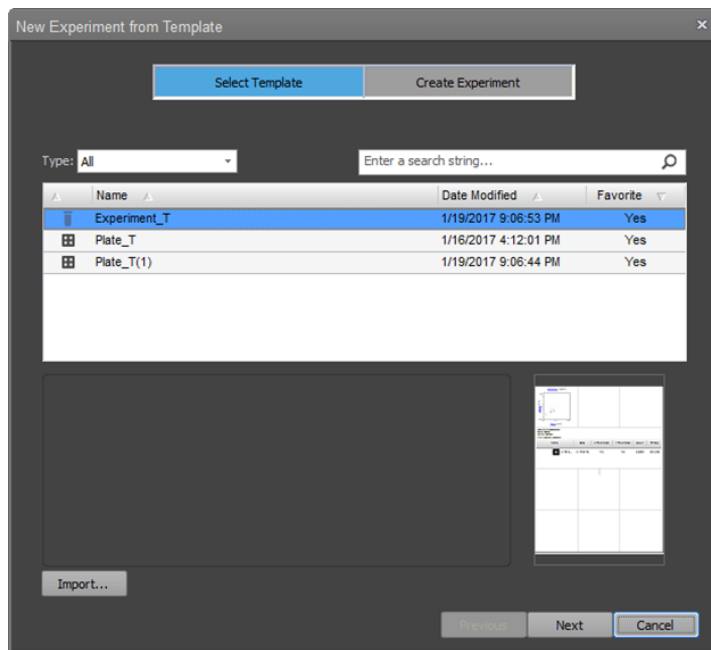


Figure 127 New Experiment from Template – Select Template

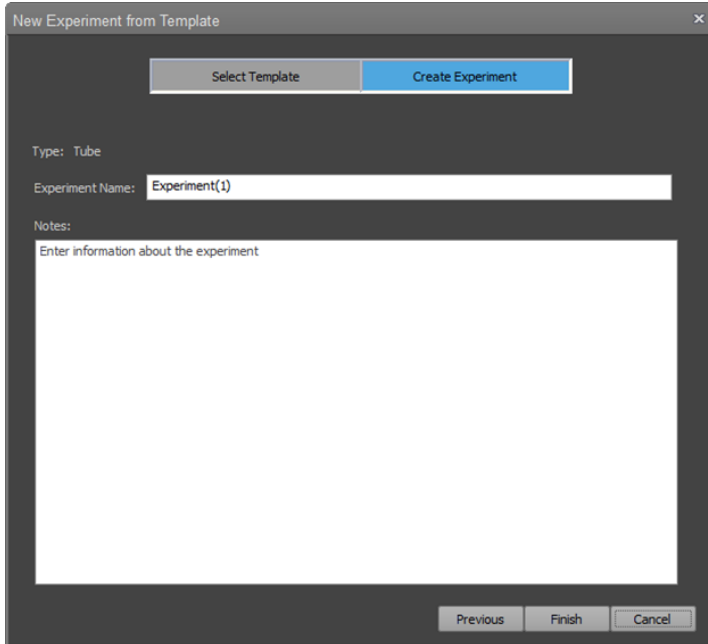
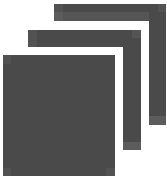


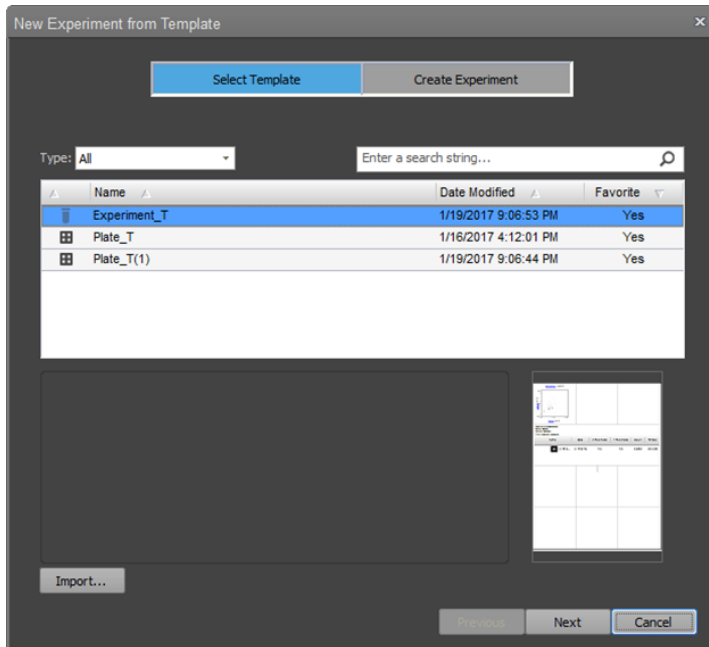
Figure 128 New Experiment from Template – Create Experiment

## Launch the new experiment from template dialog

On the Ribbon bar, click **File** to open the File menu, then select **New Experiment from Template**.



Alternatively, click the **New Experiment from Template** icon in the *Home* tab.



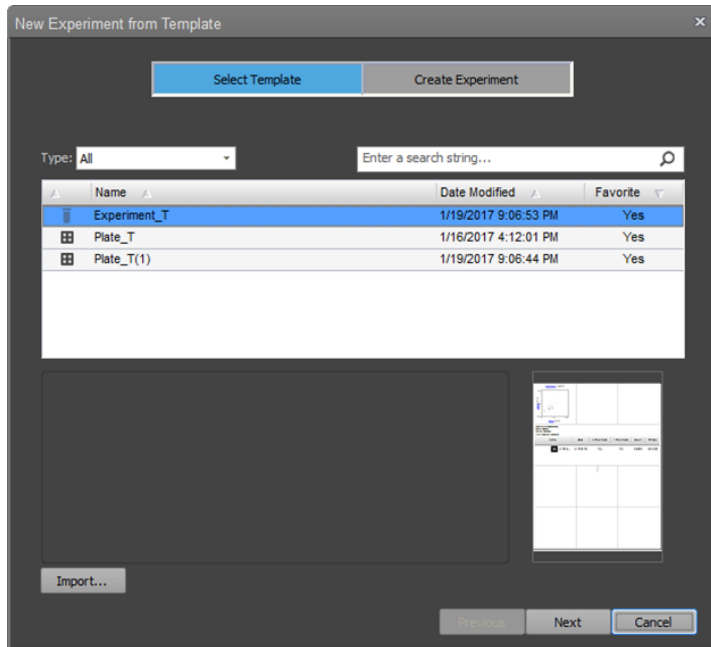
## Select template

The first step in creating a New Experiment from Template is template selection. In a template, the following Experiment settings are already defined: Workspace, Instrument settings, Run Protocols, Heat Map settings, Compensation, Overlays, Hierarchy, and the number of mapped Groups and Samples.

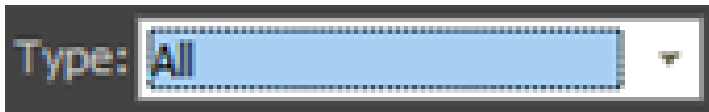
1. If it is not already selected, click **Select Template**.



Select Template tab of the New Experiment from Template dialog opens. The Select Template tab contains a list of available templates and allows you to filter the list by template type or by user-defined keywords found in the template name or description.



- To filter the template list by template type, select the desired type from the **Type** dropdown. Available options are **All**, **Tube**, or **Plate**.



The template list displays only the type of template selected in the dropdown.

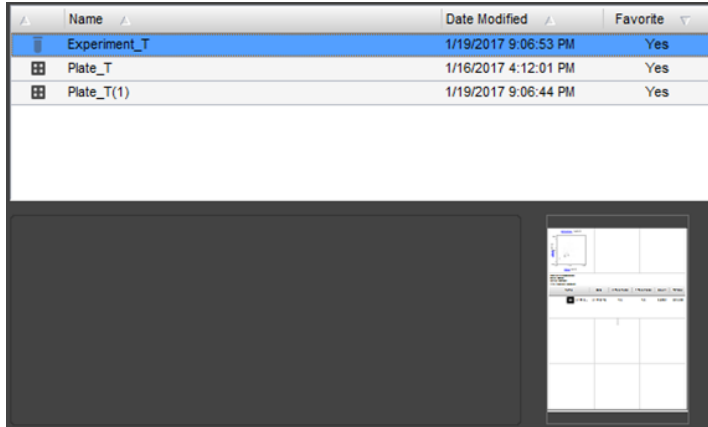
- To search the template list by keywords found in the template name or description, type in the keyword in the search text box and press **Enter**.



The template displays only the templates that contain the user-defined keyword in their name or description.

- To sort the template listed by Type, Name, or Date Modified, click on the corresponding column heading.
- To import a template that is not displayed in the template list, click **Import** to open the *File Open (Import) dialog* with the filter set to Attune™ Experiment Template (\*.aet). Select the \*.aet file that corresponds to the template you wish to add to the template list, then click **Open**. The imported template is displayed and selected in the template list.

- To select a template from the template list, click the desired template in the list. The selected template will be highlighted in the list, the description field displays the template description (if available), and the thumbnail shows the thumbnail image that corresponds to your selected template.



- To select the template for your new experiment, select the desired template in the list, then click **Next**.

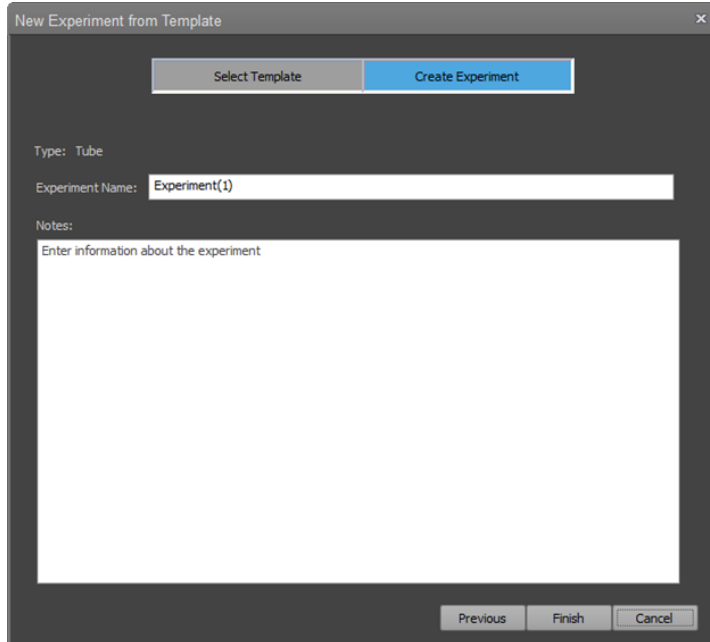
Alternatively, click the **Create Experiment** tab after selecting the template from the list.



Create Experiment tab of the New Experiment from Template dialog opens.

## Create experiment

The second step in creating a New Experiment from Template is the addition of Experiment name and Experiment notes in the Create Experiment tab.



- The Create Experiment tab displays the Experiment type selected in the Select Template tab (in this example, Tube experiment).
- By default, the Experiment Name is populated with Experiment or Plate (for Tube and Plate experiments, respectively).



1. To change the Experiment Name, click on the Experiment Name text box and type in the desired name.
  - You can enter up to 50 characters; the following characters are not permitted: \ / : \* ? " < > | %.
  - The software automatically removes trailing or leading spaces.
  - Names cannot end with a period character or be any of the following system reserved names: CON, PRN, AUX, CLOCK\$, NUL, COM1, COM2, COM3, COM4, COM5, COM6, COM7, COM8, COM9, LPT1, LPT2, LPT3, LPT4, LPT5, LPT6, LPT7, LPT8, LPT9.
  - Duplicate names are not permitted. If the name is already used, the name is appended with "(#)", where # is the next available unique number. This number increments as necessary to assure that the name is unique. For example, Experiment(2), Experiment(3), Experiment(4), etc.



2. *Optional:* Enter notes for the Experiment. You can enter up to 500 characters; any character is permitted.



3. Click **Finish** to create the New Experiment and close the dialog.
  - For Tube-only Experiments, the software opens the first Sample in the Experiment and displays the Experiment Workspace (“Workspace levels” on page 118).
  - For Plate Experiments, the software displays the Heat Map view (Chapter 6, “Heat Map View”).

To go to the Select Template tab to select another template, click **Previous**.

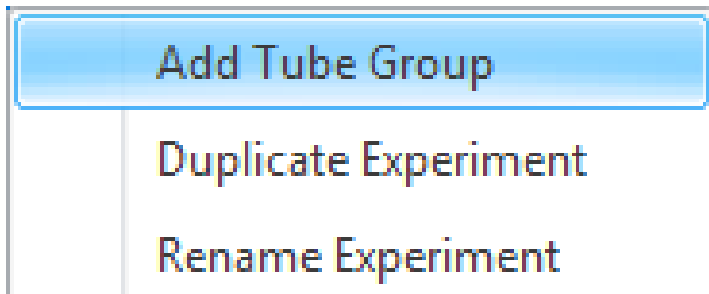
To close the dialog without creating an Experiment, click **Cancel**.

## New group dialog

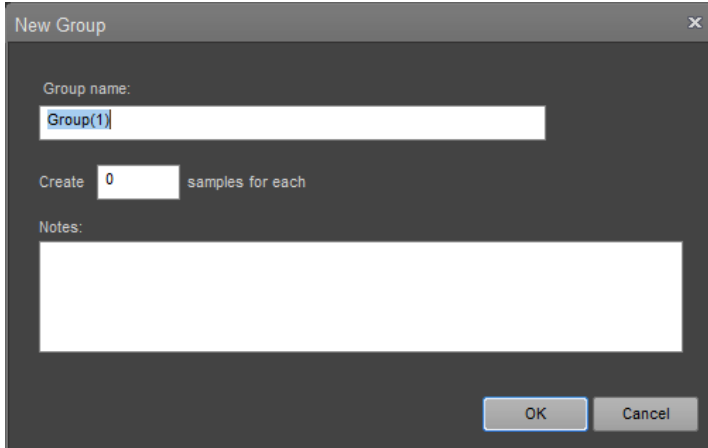
The *New Group* dialog allows you to create a new Tube group and specify the number of Samples to create for that Group.

### Launch the new group dialog

On the *Experiment Explorer*, right-click on an Experiment and select **Add Tube Group**.



*New Group* dialog opens.



## Create a group

1. Accept the default Group name or enter a new name.



- You can enter up to 50 characters; the following characters are not permitted: \ / : \* ? " < > | %.
- The software automatically removes trailing or leading spaces.
- Names cannot end with a period character or be any of the following system reserved names: CON, PRN, AUX, CLOCK\$, NUL, COM1, COM2, COM3, COM4, COM5, COM6, COM7, COM8, COM9, LPT1, LPT2, LPT3, LPT4, LPT5, LPT6, LPT7, LPT8, LPT9.
- Duplicate names are not permitted.
- The default name for a Group is **Group**. If a Group using the default name already exists, a numerical suffix is added in parentheses to ensure a unique name; for example, **Group (2)**.

2. Enter the number of Tube samples to create for the Group.



- For a Tube-only Experiment, you can enter a number from **1** to **400**. If the number of Groups times the number of Samples is  $>400$ , the number of Samples within a Group will revert to a number that is  $\leq 400$ .
  - For a 96-well plate, you can enter a number from **1** to **304**. If the number of Groups times the number of Samples is  $>304$ , the number of Samples within a Group will revert to a number that is  $\leq 304$ .
  - For a 384-well plate, you can enter a number from **1** to **16**. If the number of Groups times the number of Samples is  $>16$ , the number of Samples within a Group will revert to a number that is  $\leq 16$ .
  - The software adds the number of Samples to Experiment Explorer.
3. The software creates a unique name for each Sample with the default Sample name set in the Options dialog ► General tab and appends it with a numerical suffix; for example, **Sample (2)**.
  4. *Optional:* Enter notes for the Experiment. You can enter up to 500 characters; any character is permitted.



5. Click **OK** to create the new Group and close the *New Group* dialog.  
To close the dialog without creating a Group, click **Cancel**.

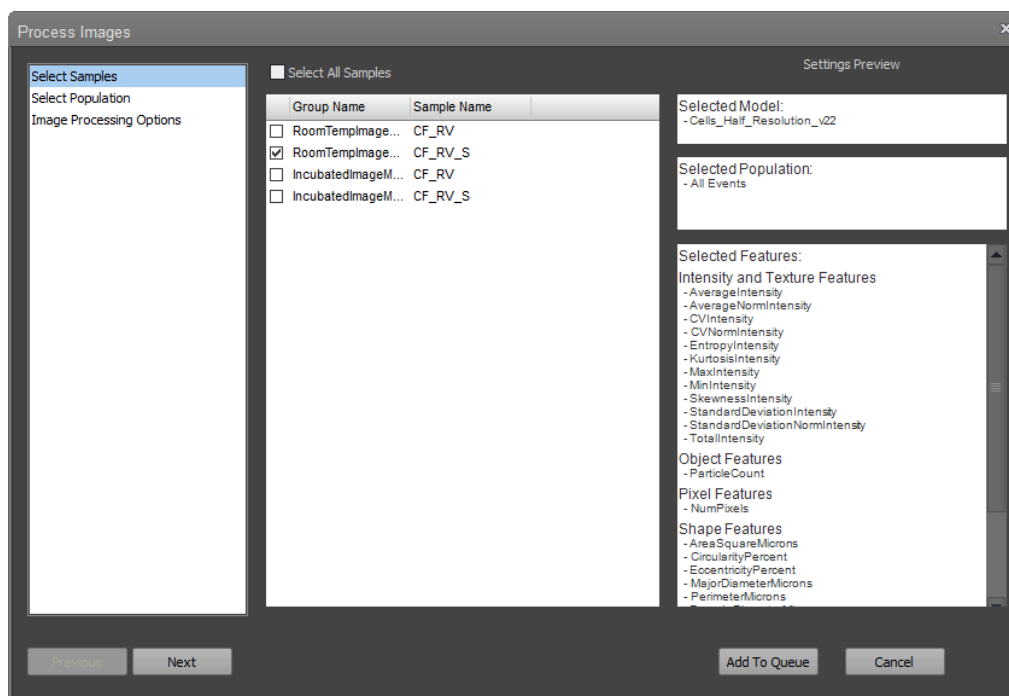
## Overview

**Process Images** dialog enables you to process captured images from selected **Samples** and **Populations** using system-provided or imported **Image Processing Models**.

---

**IMPORTANT!** Images can only be processed for Samples in the active Attune™ CytPix™ Experiment.

---



**Process Images** dialog consists of three screens:

- **Select Samples:** Enables you to select **Samples** for which to process images (see “Select Samples” on page 624).
- **Select Population:** Enables you to select one or more **populations (gates)**, where only events with the selected gates are processed (see “Select Population” on page 626).
- **Image Processing Options:** Provides options to select an **image processing model** (see “Image Processing Options” on page 628).

## Open the Process Images dialog

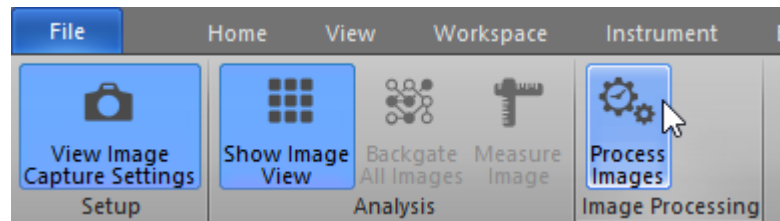
There are two ways to open the **Process Images** dialog to start image processing:

1. On the **Image Settings** ribbon tab **Image Processing** controls, click **Process Images**.

---

**Note:** The **Image Processing** controls are visible only for Attune™ CytPix™ experiments.

---

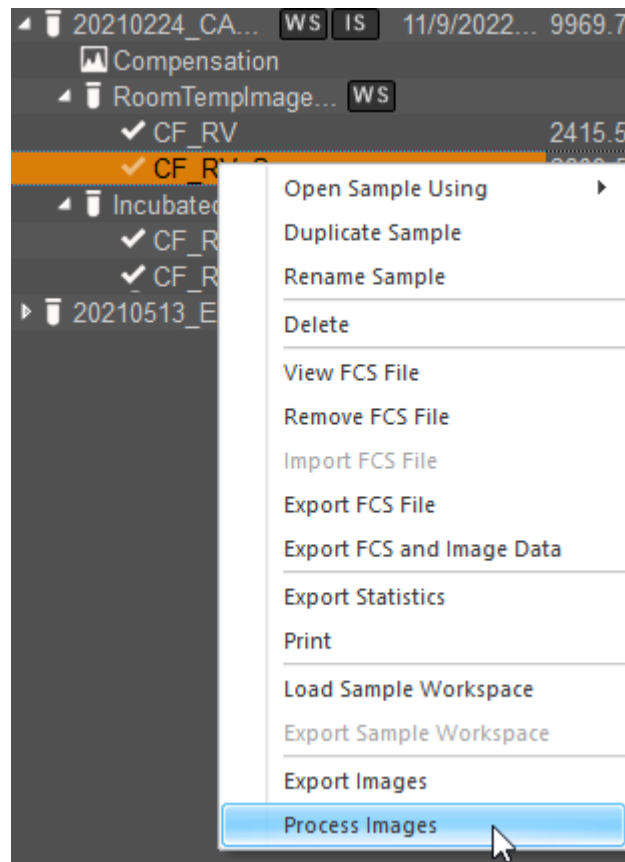


2. Alternatively, select **Process Images** in an **Experiment Explorer** context menu.

---

**Note:** **Process Images** option is only available for experiments that have image data.

---



## Select Samples

The **Select Samples** screen enables you to select the **Samples** for which to process images.

**Note:** The **Select Samples** screen is the only required screen of the **Process Images** dialog. You **must** select Samples to process images.

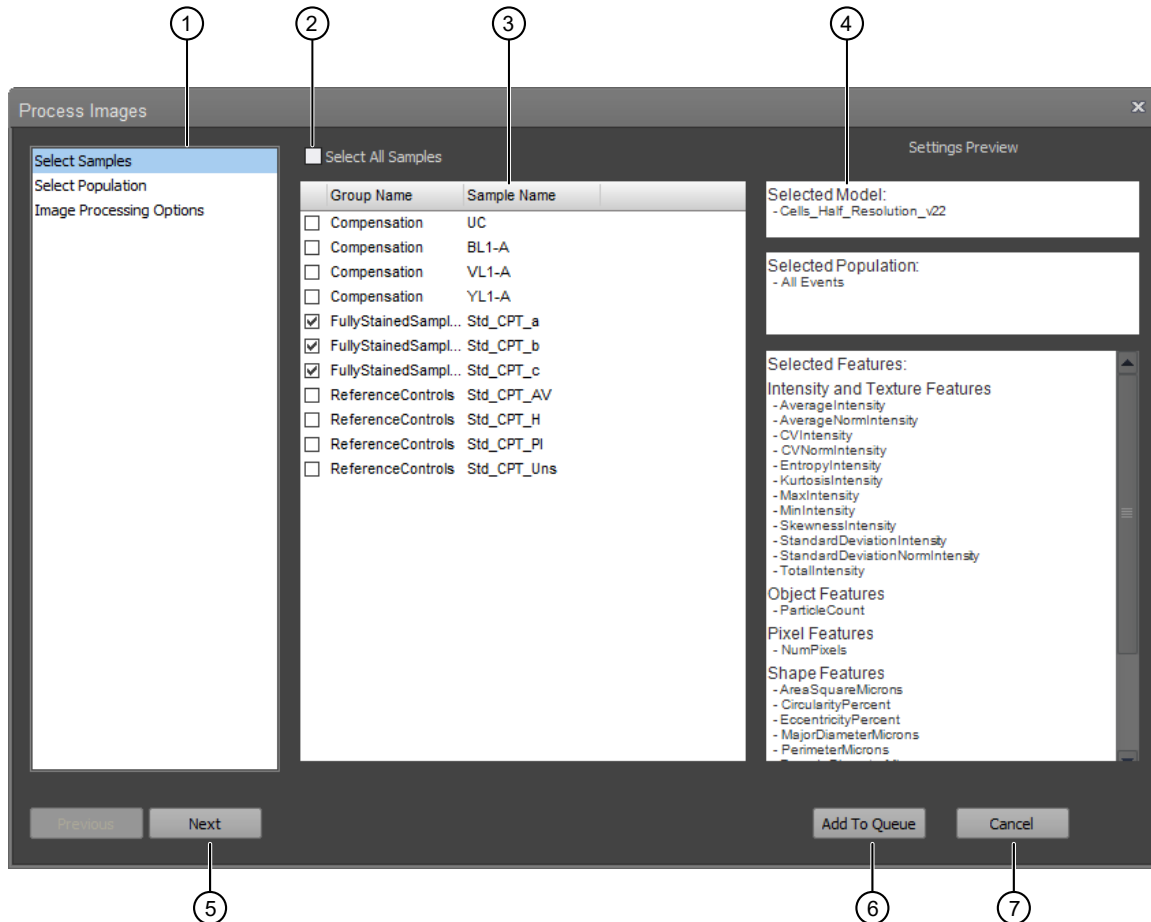
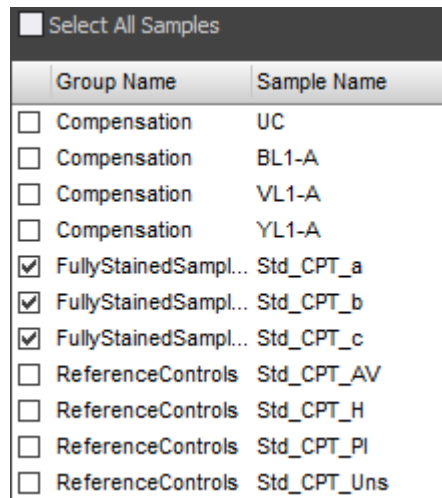


Figure 129 Select Samples screen of Process Images dialog.

- |                           |                |
|---------------------------|----------------|
| ① Dialog navigation panel | ⑤ Next         |
| ② Select All Samples      | ⑥ Add To Queue |
| ③ Sample selection list   | ⑦ Cancel       |
| ④ Settings Preview        |                |

- **Dialog navigation panel:** Selects the dialog screen that is displayed, and enables you to go between dialog screens non-sequentially.
- **Sample selection list:** Shows the available Samples in the Experiment that can be selected for image processing.
  - To select a Sample for image processing, select the corresponding **Sample checkbox**. You can select multiple Samples for processing.
  - To remove a Sample from the image processing workflow, deselect the corresponding **Sample checkbox**.

- If the **Process Images** dialog is opened from the **Experiment Explorer**, the Samples are automatically selected based on the **Sample, Group, or Experiment** that was used to open the dialog.



- To select or deselect all Samples in the Experiment, select or deselect the **Select All Samples** option above the sample selection list.
- **Next:** Opens the next screen of the dialog (**Select Population** screen)
- **Add To Queue:** Adds the Samples to the image processing queue to be processed serially in the order they were added (see “Image processing queue” on page 630).
- **Cancel:** Closes the dialog without adding the Samples to the processing queue.

## Select Population

The **Select Population** screen enables you to select one or more populations (gates) where only events within the gates are processed. The data shown and list of gates is based on the active workspace when the **Process Images** dialog is opened.

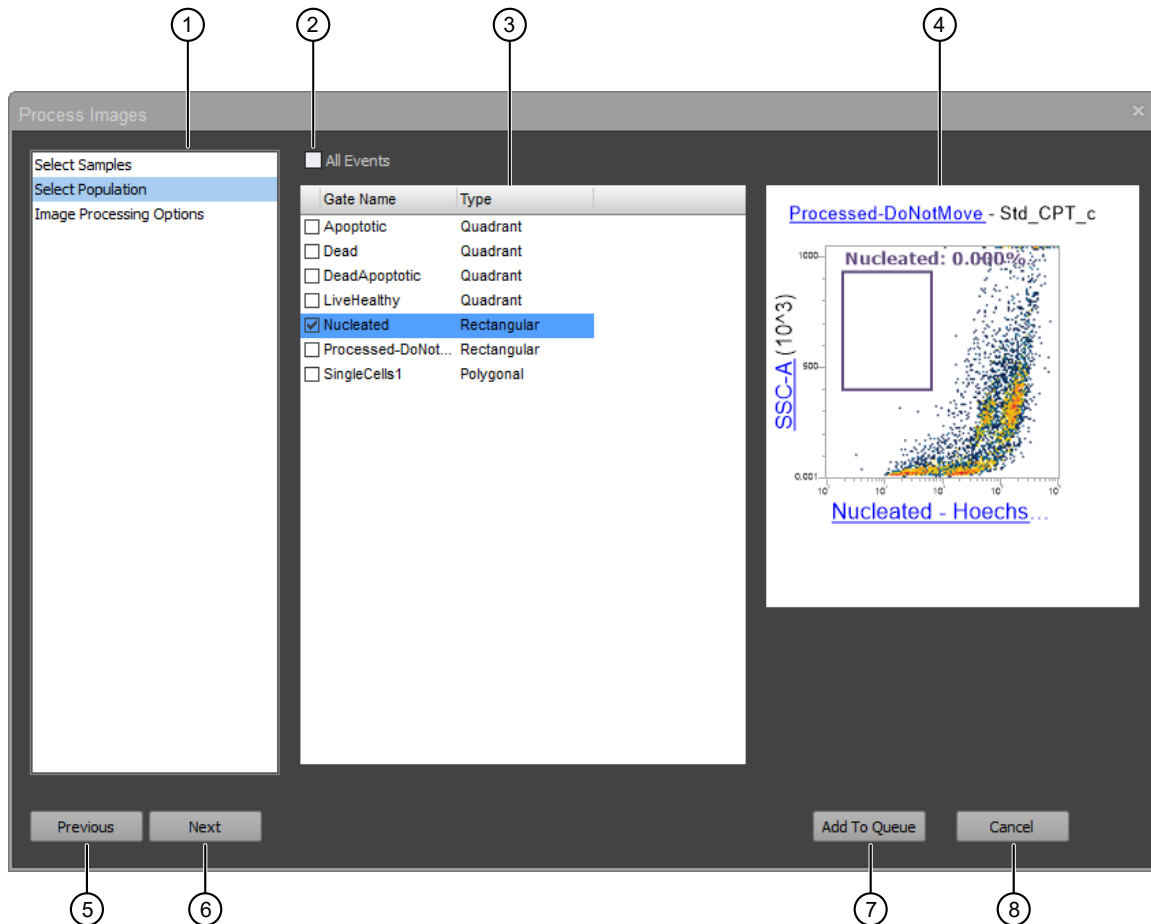


Figure 130 Select Population screen of Process Images dialog

- |                             |                |
|-----------------------------|----------------|
| ① Dialog navigation panel   | ⑤ Previous     |
| ② All Events                | ⑥ Next         |
| ③ Population selection list | ⑦ Add To Queue |
| ④ Preview plot              | ⑧ Cancel       |

- **Dialog navigation panel:** Selects the dialog screen that is displayed, and enables you to go between dialog screens non-sequentially.
- **Population selection list:** Shows the available populations (gates) that can be selected to filter the images to be processed. By default, **All Events** is selected.
  - To select a population for image processing, select the corresponding **Population checkbox**. You can select multiple populations.
  - To remove a population from the image processing workflow, deselect the corresponding **Sample checkbox**.



- To select or deselect all events in the Sample for image processing, select or deselect the **All Events** option above the population selection list. By default, **All Events** is selected.
- The Attune™ Cytometric Software uses the selected gate to process all Samples. If the gate does not exist on the selected Workspace, the software tries the next level up on the gating hierarchy (Sample>Group>Experiment), until it cannot find the gate and reverts to **All Events** for that Sample.
- **Preview plot:** Shows the selected gate and FCS data of the active sample. When no gate is selected, the preview plot does not show any plot.
- **Previous:** Opens the previous screen of the dialog (**Select Sample** screen).
- **Next:** Opens the next screen of the dialog (**Select Population** screen).
- **Add To Queue:** Adds the Samples to the image processing queue to be processed serially in the order they were added (see “Image processing queue” on page 630).
- **Cancel:** Closes the dialog without adding the Samples to the processing queue.

## Image Processing Options

The **Image Processing Options** screen enables you to select the **Image Processing Model** to use and set the selected model as the default. It also lists the **Image Processing Parameters** for the selected model.

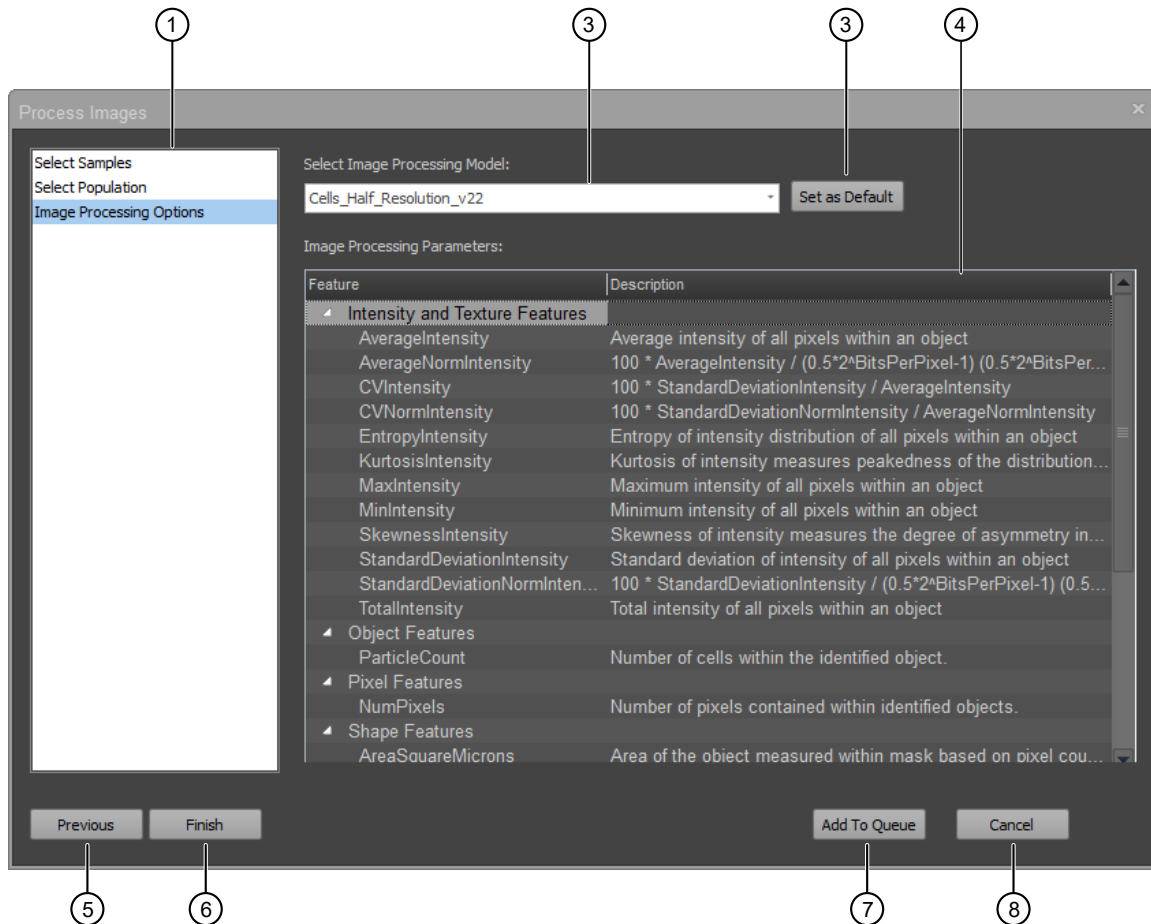


Figure 131 Image Processing Options screen of Process Images dialog

- |                                 |                |
|---------------------------------|----------------|
| ① Dialog navigation panel       | ⑤ Previous     |
| ② Select Image Processing Model | ⑥ Finish       |
| ③ Set as Default                | ⑦ Add To Queue |
| ④ Image Processing Parameters   | ⑧ Cancel       |

- **Dialog navigation panel:** Selects the dialog screen that is displayed, and enables you to go between dialog screens non-sequentially.
- **Select Image Processing Model:** Enables you to select an **Image Processing Model** to use. By default, **All Events** is selected.
  - Click the **Select Image Processing Model** dropdown, then select the model you want to use to process the selected images. Currently available image processing models are: **Beads Only (Full Resolution)**, **Cells (Full Resolution)**, **Cells (Half Resolution)**.



- Click **Set As Default** to set the selected image processing model as the default for subsequent image processing operations.

---

**Note:** You can also change the default image processing model in the **Options dialog ▶ Image Options tab** using the **Set as favorite** button (see “Image Options” on page 673).

- The Attune™ Cytometric Software uses the selected image processing model to process the selected Samples.

---

**IMPORTANT!** Regardless of the size of the captured images in the experiment, images are processed at either full resolution (248 × 248 pixels) or at half resolution (124 × 124 pixels). Processing the images at half resolution decreases data footprint by 4-fold and improves the processing times by >4-fold. However, this comes with a potential trade off in accuracy.

- **Image Processing Parameters:** Shows the list of the parameters that are generated when the image processing is completed and provides a description of each parameter.

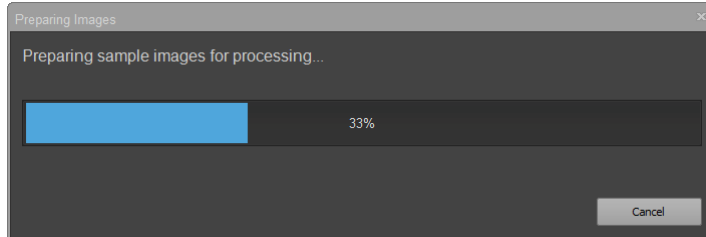
---

**Note:** For a list and description of image processing parameters, see Appendix B, “Attune™ Cytometric Software image processing parameters”. The Attune™ Cytometric Software does not allow user-selected image processing parameters.

- **Previous:** Opens the previous screen of the dialog (**Select Population** screen).
- **Finish:** Adds the Samples to the image processing queue to be processed serially in the order they were added (see “Image processing queue” on page 630).
- **Add To Queue:** Adds the Samples to the image processing queue to be processed serially in the order they were added.
- **Cancel:** Closes the dialog without adding the Samples to the processing queue.

## Image processing queue

- When **Add to Queue** or **Finish** is clicked on the **Process Images** dialog, all selected samples are added to a processing queue and are processed serially in the order they were added.
- The software displays the progress of the image processing operation as Samples are added to the queue and gated events are calculated.



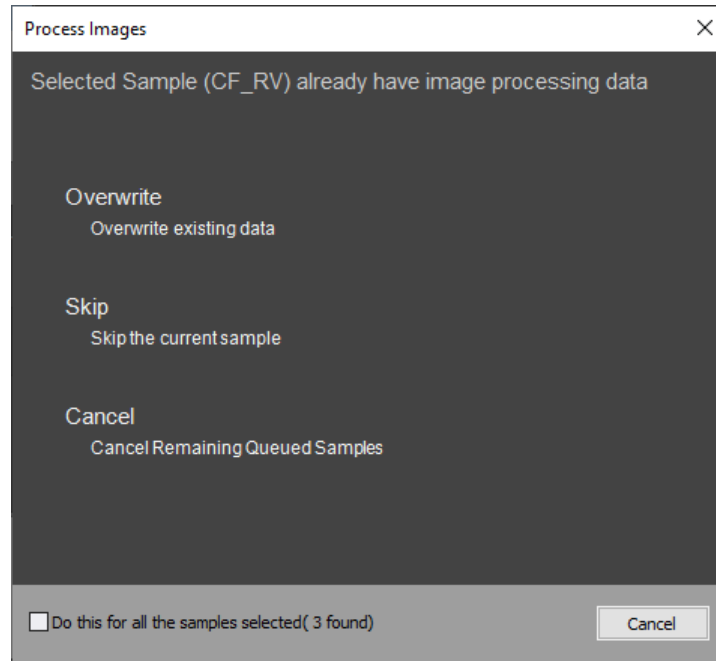
- You can manage the sample queue using the Attune™ Image Processing Dashboard (see Chapter 24, “Attune™ Image Processing Dashboard”).

Attune Image Processing Dashboard

Manage Queue		Search						
Scheduling	Batch ID	Experiment Name	Group Name	Sample Name	Model	Status	Estimated Time	Progress
Configuration	2022-11-17-1 2-6-3	20210224_CARTRam os	IncubatedImage More	CF_RV	Cells_Half_Resolution_v22	In Progress	17.3 Min	9 %
	2022-11-17-1 2-6-54	20210224_CARTRam os	IncubatedImage More	CF_RV_S	Cells_Half_Resolution_v22	Pending	0.0 Min	0 %

- Image processing proceeds even if you close the Attune™ Cytometric Software.
- Image processing is paused during sample acquisition.
- If a specific population was selected in the **Select Population** screen, the events contained in each selected gate are calculated for each Sample.  
If no gates were selected in the **Select Population** screen, the software reverts to **All Events** for the Sample and processes all captured images in the Sample.
- If no image processing model is selected in the **Image Processing Options** screen, the software uses the default image processing model.
- You can set an image processing model as the default for subsequent image processing operations by selecting it in the **Image Processing Options** screen, then clicking **Set As Default** (see “Image Processing Options” on page 628).  
Alternatively, you can select the desired image processing model in the **Options dialog** ▶ **Image Options tab**, then click **Set as favorite**. (see “Image Options” on page 673).

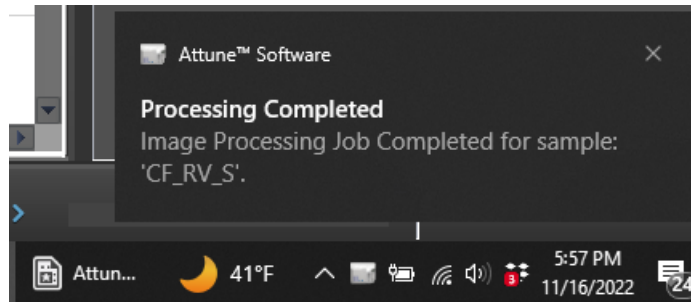
- If the selected samples already have image processing data, the software displays a warning dialog that enables you to overwrite the existing data, to skip the current sample, or to cancel image processing (see “Reprocess images” on page 634).



## Completion of image processing

When the Attune™ Cytometric Software is running and image processing is completed for any sample in the processing queue that is specific to the currently logged-in user, the software provides the following notifications:

- System tray displays the **Processing Complete** notification.

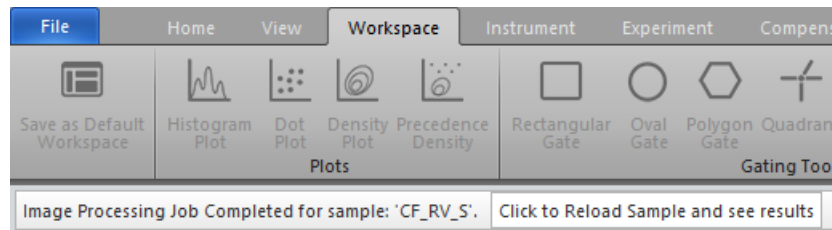



---

**Note:** Windows notifications must be enabled to see system tray notifications.

---

- Message bar displays the **Image Processing Job Complete** notification with an option to reload the sample and see the results.




---

**Note:** Clearing the message by clicking **X** or **Reload Sample and see results** button hides the message bar.

---

- If the Attune™ Software is running in the SAE ("Security, Auditing, and Electronic Signature") mode and the experiment was signed, the message bar displays a warning that indicates that changes to the experiment will invalidate any existing signatures.

When image processing is completed for a sample, the icon next to the sample in the Experiment Explorer is updated to indicate that the sample has both FCS and image processing data.

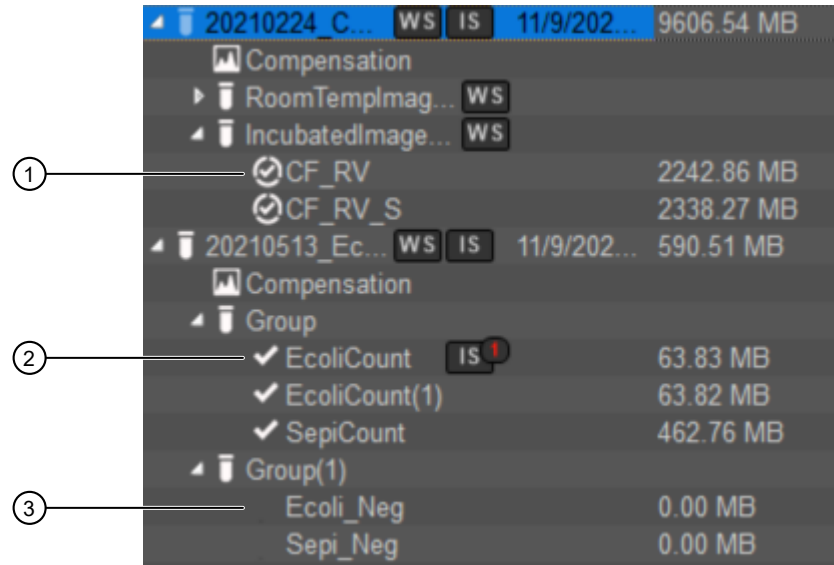


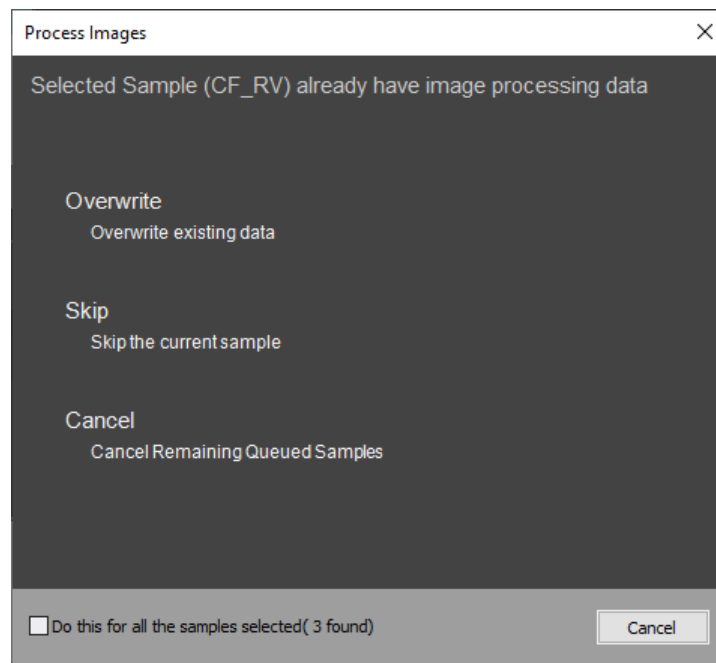
Figure 132 Experiment Explorer containing samples with and without image processing data

- ① Sample with FCS and Image Processing data
- ② Sample with FCS data only
- ③ Sample with no data

## Reprocess images

You can reprocess images in a sample to allow a different processing model or algorithm, or to use other populations or features to extract cell image information. When samples are reprocessed, any ongoing processing is canceled and the new processing request is added to the queue.

1. Select the samples, populations, and image processing options as described in Process images. When you start processing a request for a sample that has already been processed, the software displays a dialog that states that the selected sample already has image processing data. The software displays the dialog for each sample that meets this condition.



2. To process the selected samples and overwrite the existing image processing data, select **Overwrite**.

---

**Note:** If you have selected multiple samples with existing image processing data, you can apply the **Overwrite** option to all samples that meet this condition by selecting **Do this for all the samples selected** checkbox. When the last sample that meets this condition is confirmed, the image processing dialog is closed.

---

The selected samples that already have image processing data are added to the queue and processed along with any samples that do not have existing data.

---

**IMPORTANT!** When image processing completes for each sample, any existing results are overwritten.

---



3. To skip the samples with existing image processing data and process only the samples that do not have existing data, select **Skip**.

---

**Note:** If you have selected multiple samples with existing image processing data, you can apply the **Skip** option to all samples that meet this condition by selecting **Do this for all the samples selected** checkbox. When the last sample that meets this condition is confirmed, the image processing dialog is closed.

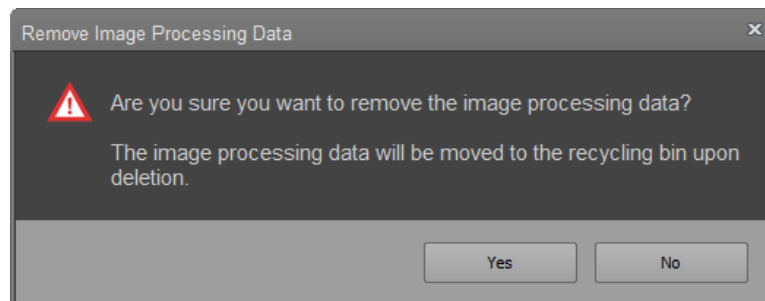
---

The selected samples that already have image processing data are not added to the queue and only those samples that do not have existing data are processed.

4. To cancel the processing request without adding the selected samples to the image processing queue, select **Cancel**.  
The image processing dialog remains open, allowing you to update the sample selection, if desired.

## Remove Image Processing Data dialog

The **Remove Image Processing Data dialog** helps to prevent image processing data from being accidentally removed from Samples. The dialog is displayed when you select **Remove Image Processing Data** option from a **Sample context menu**.



- To remove the image processing data from the sample, click **Yes**. The image processing data are moved to the recycling bin.
- To close the dialog without removing the image processing data, click **No** or **X**.

## Attune™ Image Processing Dashboard

The Attune™ Image Processing Dashboard is automatically opened when you power on the computer that runs the Attune™ Cytometric Software and it functions independently of the Attune™ Software.

The Attune™ Image Processing Dashboard enables you to:

- View image processing queue, status, and progress (page 638)
- Reprioritize the image processing queue (page 638)
- Cancel image processing requests (page 638)
- Pause or resume image processing (page 638)
- Schedule image processing



Figure 133 Attune™ Image Processing Dashboard

- ① Manage Queue tab
- ② Image processing queue
- ③ Move request up or down buttons
- ④ Delete button
- ⑤ Resume processing button
- ⑥ Pause processing button
- ⑦ Close dashboard button

## View image processing queue, status, and progress

You can view the image processing queue, status, and progress in the **Manage Queue** tab of the Attune™ Image Processing Dashboard:

	Batch ID	Experiment Name	Group Name	Sample Name	Model	Status	Estimated Time	Progress
Scheduling	2022-11-17-1	20210224_CARTRam	IncubatedImage	CF_RV	Cells_Half_Resolution_v22	In Progress	17.3 Min	9 %
Configuration	2-6-3	os	More					
	2022-11-17-1	20210224_CARTRam	IncubatedImage	CF_RV_S	Cells_Half_Resolution_v22	Pending	0.0 Min	0 %
	2-6-54	os	More					

Figure 134 Image processing queue

- **Batch ID:** Unique identification assigned to the image processing job based on the date and time the sample was added to the processing queue.
- **Experiment Name:** Name of the experiment in Experiment Explorer from which the sample was selected for image processing.
- **Group Name:** Group name in Experiment Explorer from which the sample was selected for image processing.
- **Sample Name:** Name of the sample in Experiment Explorer that was selected for image processing.
- **Model:** Image processing model selected in Process Images wizard.
- **Status:** Image processing status for the sample (In Progress, Pending, or Paused).
- **Estimated Time:** Estimated time for the completion of image processing for the sample.
- **Progress:** Percentage of the image processing job that is completed.

---

**Note:** Image processing is paused during sample acquisition.

---

## Reprioritize the image processing queue

To move a processing request up or down in the image processing queue, select the request, then use the **Move to Top**, **Move Up**, **Move Down**, and **Move to Bottom** buttons to move it to the desired position in the queue.

## Cancel image processing requests

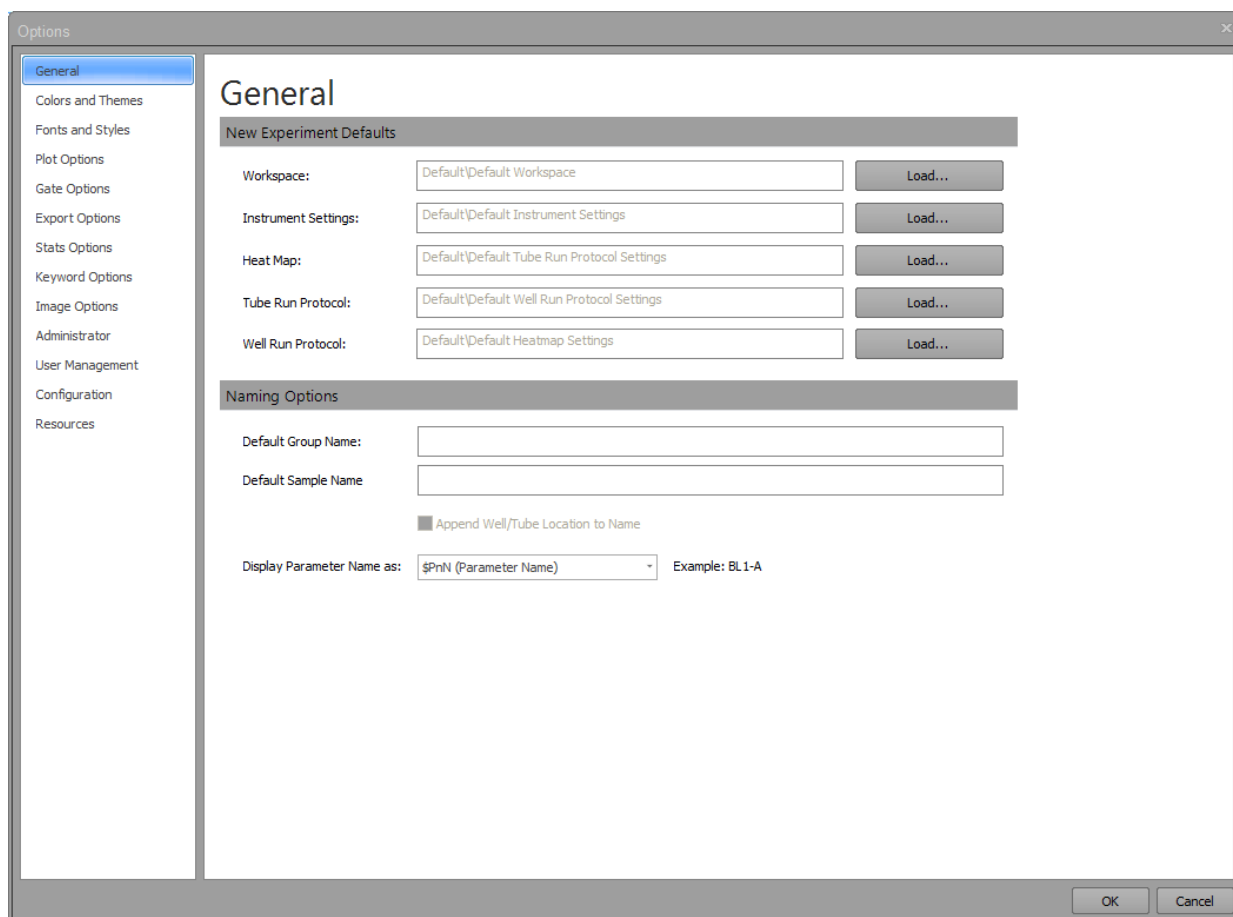
To cancel an image processing request, select the request in the image processing queue, then click **Delete**.

## Pause or resume image processing

1. To pause an ongoing image processing of a sample, select the sample, then click **Pause**.
2. To resume image processing, click **Resume**.

## Overview

The **Options dialogs** let you to customize the Attune™ Cytometric Software by configuring personal settings and changing the default options.



- To open the **Options dialog**, click the **Options** button on the **Quick Access toolbar** (“Quick Access toolbar” on page 52). This option is enabled only when the instrument is not acquiring.
- The **Options dialog** is divided into tabs, which are listed on the left side of the window. Select a tab to open the tab content on the right side of the window.
- Some options are user-specific, while others are application-specific (i.e., global to all users) and customizable only by an authorized user.

For example, the **Plate Options** tab (“Plate Options” on page 675) is available only when a CytKick™ Max™ Autosampler is connected to the Attune™ instrument. Similarly, the **SAE Configuration** controls (“SAE Configuration” on page 680) in the **Administrator** tab are only visible if the SAE-specific DESKey device is present.

- The **Options dialog** contains the following tabs.
  - **General** (“General” on page 641)
  - **Colors and Themes** (“Colors and Themes” on page 647)
  - **Fonts and Styles** (“Fonts and Styles” on page 649)
  - **Plot Options** (“Plot Options” on page 653)
  - **Gate Options** (“Gate Options” on page 656)
  - **Export Options** (“Export Options” on page 662)
  - **Stats Options** (“Stats Options” on page 665)
  - **Keyword Options** (“Keyword Options” on page 669)
  - **Image Options** (“Image Options” on page 673)
  - **Plate Options** (“Plate Options” on page 675)
  - **Administrator** (“Administrator” on page 676)
  - **User Management** (“User Management” on page 687)
  - **Configuration** (“Configuration” on page 695)
  - **Resources** (“Resources” on page 701)
- By default, the **General** tab is displayed when the **Options dialog** is first opened.
- When you make changes in the options menu, click **OK** to save the changes and close the **Options dialog**.

Click **Cancel** or the **X** in the top right corner to close the dialog and revert to the previous selections without applying the changes.
- When you switch between tabs or click **OK**, all fields are validated.

Any invalid fields in the currently selected tab view are indicated on a field-by-field basis. You cannot switch views or complete the **OK** action until all errors are corrected.

# General

## Overview

**General** tab is selected by default when the **Options dialog** is first opened. It enables you to view the **New Experiment Defaults** and to change the **Naming Options** for Groups, Samples, and Display Parameters (“Naming Options” on page 644).

The screenshot shows the 'Options' dialog box with the 'General' tab selected. The dialog is divided into two main sections: 'New Experiment Defaults' and 'Naming Options'.

**New Experiment Defaults:**

- Workspace: Default\Default Workspace (with a 'Load...' button)
- Instrument Settings: Default\Default Instrument Settings (with a 'Load...' button)
- Heat Map: Default\Default Tube Run Protocol Settings (with a 'Load...' button)
- Tube Run Protocol: Default\Default Well Run Protocol Settings (with a 'Load...' button)
- Well Run Protocol: Default\Default Heatmap Settings (with a 'Load...' button)

**Naming Options:**

- Default Group Name: [Text input field]
- Default Sample Name: [Text input field]
- Append Well/Tube Location to Name
- Display Parameter Name as: \$PnN (Parameter Name) (dropdown menu) Example: BL1-A

At the bottom right of the dialog are 'OK' and 'Cancel' buttons.

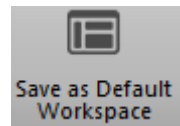
## New Experiment Defaults

**New Experiment Defaults** indicates the location of default files for **Experiment Workspace**, **Instrument Settings**, **Heat Map** settings, **Tube Run Protocol**, and **Well Run Protocol**. No changes can be made to the defaults using this menu.

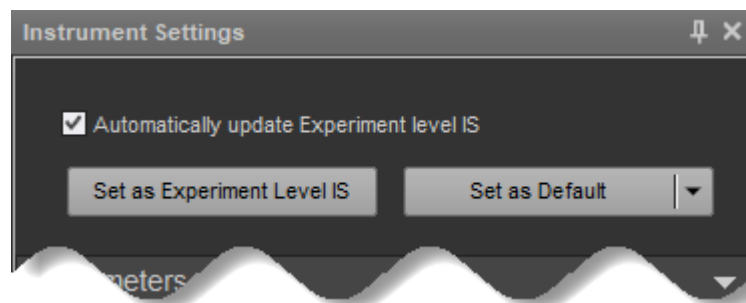
New Experiment Defaults		
Workspace:	Default\Default Workspace	Load...
Instrument Settings:	Default\Default Instrument Settings	Load...
Heat Map:	Default\Default Tube Run Protocol Settings	Load...
Tube Run Protocol:	Default\Default Well Run Protocol Settings	Load...
Well Run Protocol:	Default\Default Heatmap Settings	Load...

- **Workspace** indicates the default **Experiment-level Workspace** for each new Experiment. The **Load** button is inactive and cannot be clicked.

To change the default **Workspace**, save a **Workspace** as the default using the **Save as Default Workspace** button in the **Workspace** tab of the **Ribbon bar** (see “Save as Default Workspace” on page 81). Default Workspaces are user-account specific.



- **Instrument Settings** indicates the default **Experiment-level Instrument Settings (IS)** for each new Experiment. The **Load** button is inactive and cannot be clicked. To change the default Experiment-level instrument settings, click **Set as Default** in the **Instrument Settings panel** (Chapter 13, “Instrument settings panel”).



If the instrument configuration is changed, the default is reset to the default configuration for the user.

---

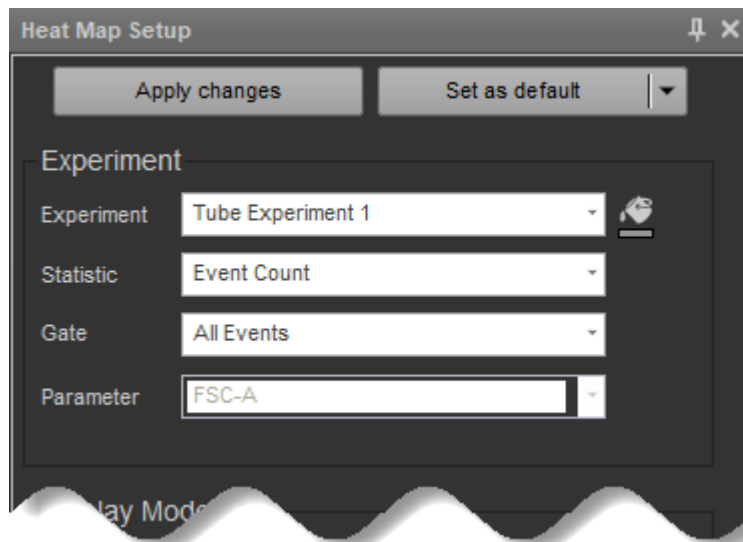
**Note:** Default **Instrument Settings (IS)** are user-account and instrument configuration specific. Changes made to the default IS of one user are not applied to the default IS of other users, except when the changes were made to system-level Instrument Settings by the Administrator.

---



- **Heat Map** indicates the default **Heat Map Settings** used for each new Experiment. The **Load** button is inactive and cannot be clicked.

To change the default **Heat Map**, click **Set as default** in the **Heat Map Setup panel** (Chapter 14, “Heat map setup panel”).

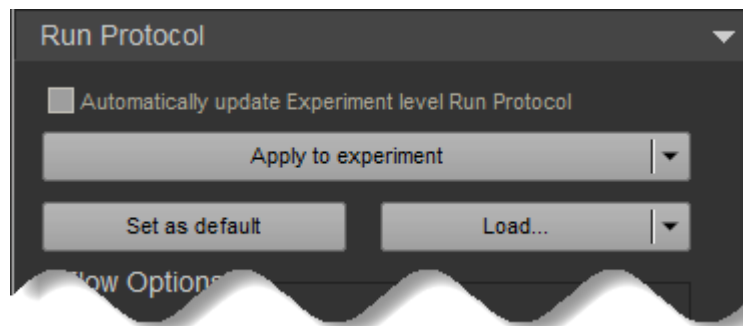


If changes are made to the default **Heat Map settings**, the default setting is changed for only that user account.

If an Experiment is created using a default Workspace that does not contain gates used to define the **Heat Map settings**, the Heat Map settings revert to **All Events** for statistics.

- **Tube Run Protocol** indicates the default **Run Protocol** to use for all new **Tube Samples** in new **Plate** or **Tube Experiments**. The **Load** button is inactive and cannot be clicked.

To change the default **Run Protocol**, click **Set as default** in the **Run Protocol** section of the **Collection Panel** (“Run protocol” on page 368).



- **Well Run Protocol** indicates the default **Run Protocol** to use for all new **Well Samples** (including **Manual Wells**) in new **Plate Experiments**. The **Load** button is inactive and cannot be clicked.

To change the default **Run Protocol** for a **Plate Experiment**, click **Set as default** in the **Run Protocol** section of the **Collection Panel** (“Run protocol” on page 368).

## Naming Options

**Naming Options** enable you to define the default Group and Sample names, and to change the way the **Display Parameter Name** is shown. Changes made to the **Default Group Name** and **Default Sample Name** are only applied to newly created Groups or Samples.

Naming Options

Default Group Name:

Default Sample Name:

Append Well/Tube Location to Name

Display Parameter Name as:  Example: BL1-A

### Default Group Name

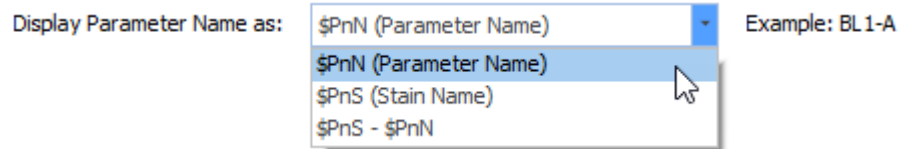
- By default, the **Default Group Name** field is blank, which sets the default Group name to **Group** when creating new Groups.
- You can enter a new name in the **Default Group Name** field, which then becomes the default Group name for all new Groups.
- Duplicate names are not permitted. If the entered name already exists, a numerical suffix in parentheses is added to the Group name; for example, **Group(2)**.
- You can enter up to 50 alpha-numeric characters in the **Default Group Name** field.
- On validation, the software removes leading and trailing spaces, and converts consecutive spaces to single spaces.
- If you attempt to enter invalid characters, a warning dialog indicates the error condition, and the invalid characters do not appear in the text field.  
The following characters are not allowed: \ / : \* < > | ?.

### Default Sample Name

- By default, the **Default Sample Name** field is blank, which sets the default Sample name to **Sample** when creating new Tube samples and to the **Well location** (e.g., **A1**) when creating new Well samples.
- You can enter a new name in the **Default Sample Name** field, which then becomes the default Sample name for all new Samples.
- Duplicate names are not permitted. If the entered name already exists, a numerical suffix in parentheses is added to the Sample name; for example, **Sample(2)**.
- You can enter up to 50 alpha-numeric characters in the **Default Sample Name** field.
- On validation, the software removes leading and trailing spaces, and converts consecutive spaces to single spaces.
- If you attempt to enter invalid characters, a warning dialog indicates the error condition, and the invalid characters do not appear in the text field.  
The following characters are not allowed: \ / : \* < > | ?.

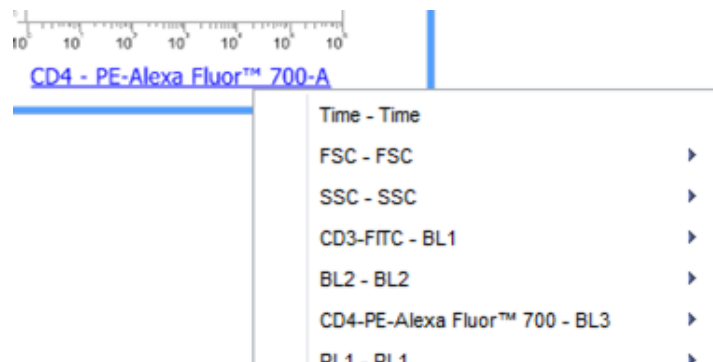
## Display Parameter Name as

- By default, the **Display Parameter Name as** is set to **\$PnN (Parameter Name)**.
- The **Display Parameter Name as** control opens a dropdown list that includes **\$PnN (Parameter Name)**, **\$PnS (Stain Name)**, and **\$PnS-\$PnN** options.

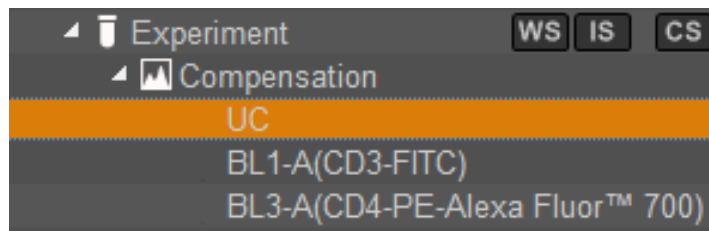


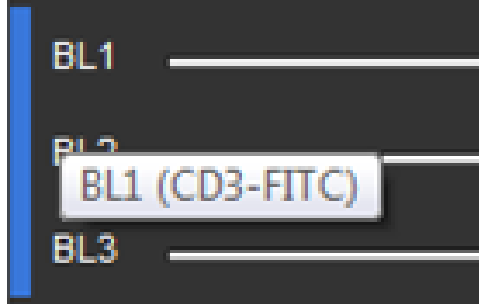
Option	Example label
\$PnN (Parameter Name)	“Example: BL1-A”
\$PnS (Stain Name)	“Example: FITC”
\$PnS-\$PnN	“Example: CD4-FITC BL1-A”

- **Display Parameter Name as** setting applies to all Experiments and affects the parameter name displayed in the **Parameter dropdown menu** for all plots, and the name displayed when the **Axis Label** checkboxes are all unchecked and no custom label has been entered.



- **Display Parameter Name as** setting also applies to the Tube or Well names for Compensation controls and to the name displayed in the tool tip for a parameter.

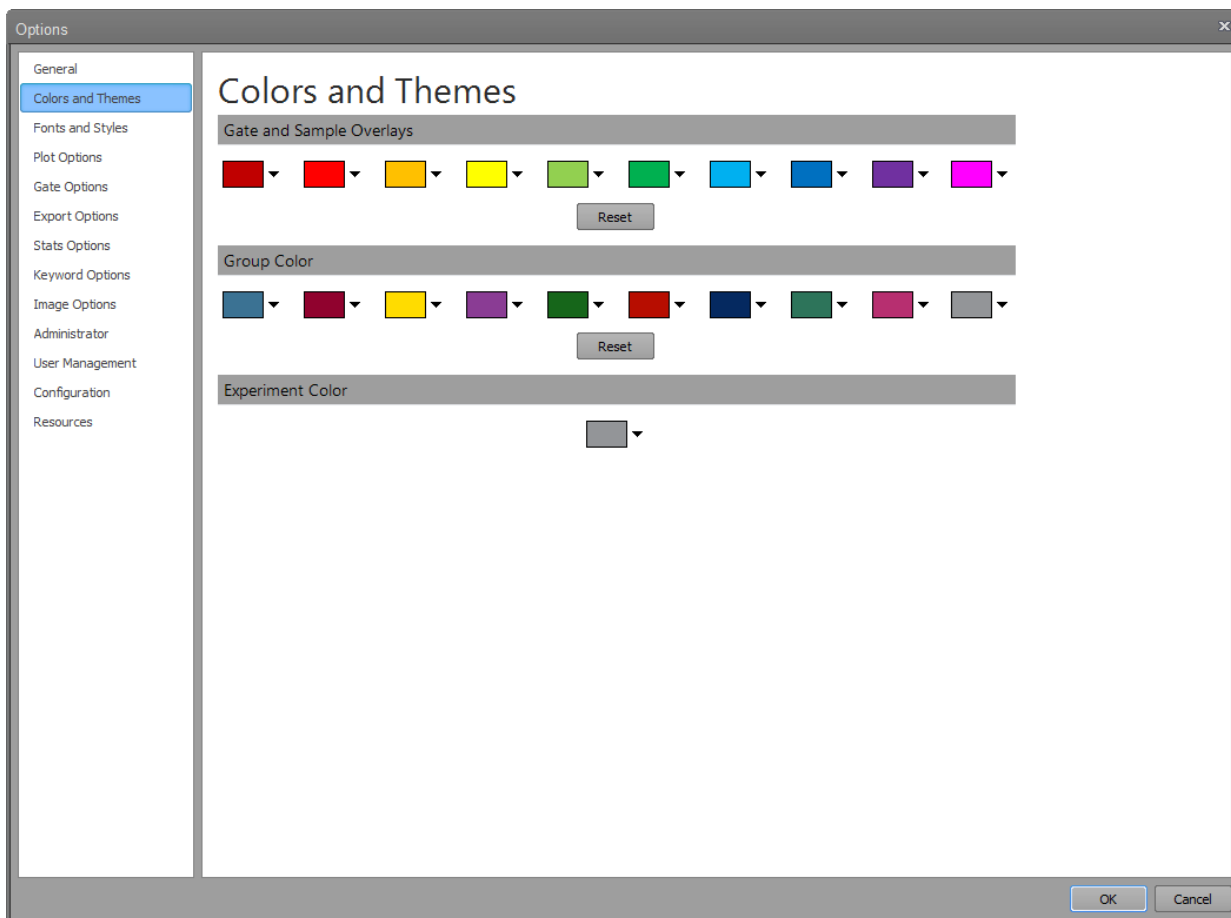




# Colors and Themes

## Overview

**Colors and Themes** lets you define the default colors for Gates, Overlays, and Experiment and Group labels.



## Colors and Themes options

- The **Colors and Themes** tab contains three groups of controls: **Gate and Sample Overlays**, **Group Color**, and **Experiment Color**.
- Gate and Sample Overlays share the defaults, but the defaults for Experiment and Group labels are controlled separately.

- Each set of defaults contains 10 standard colors. You can modify each color by clicking on the **color**, which opens the standard color picker dialog.



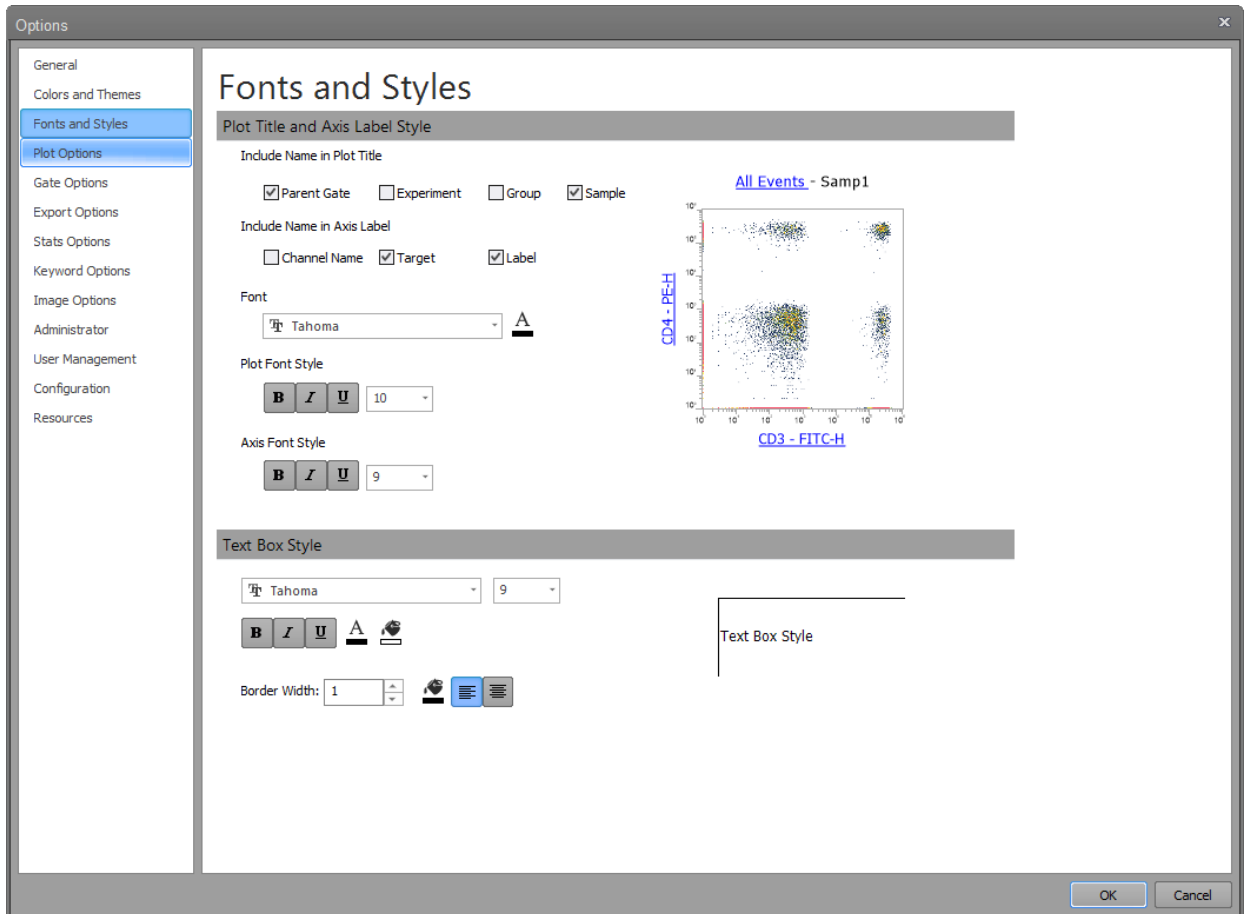
When a color is selected from the color picker dialog, the color displayed in the color dropdown is updated and the color picker dialog is closed.

- The **Experiment Color** section contains a single color dropdown with the same behavior as described above.
- Changes made to the default colors only apply to newly created Gates, Groups, and Experiments after the change is accepted.
- The **Reset** button restores the colors to the original default application colors. This option is not available on the **Experiment Color** section.

# Fonts and Styles

## Overview

**Fonts and Styles** enables you to define default display options for the text in plot labels and text boxes. The tab is divided into two sections: **Plot Title and Axis Label Style** and **Text Box Style**.



- The current settings are shown in the **Previews area** for plot and text objects on the right side of the dialog.
- Changes made in **Fonts and Styles** are applied only to newly created plots except the **Axis Label**. When **OK** is clicked, the **Axis Label** style is applied immediately and affects all past, present, and future instances.

## Plot Title and Axis Label Style

**Plot Title and Axis Label Style** controls enable you to modify the default settings for the naming format, font, and style of the plot title and axis labels.

**Plot Title and Axis Label Style**

**Include Name in Plot Title**

Parent Gate  
  Experiment  
  Group  
  Sample

**Include Name in Axis Label**

Channel Name  
  Target  
  Label

**Font**

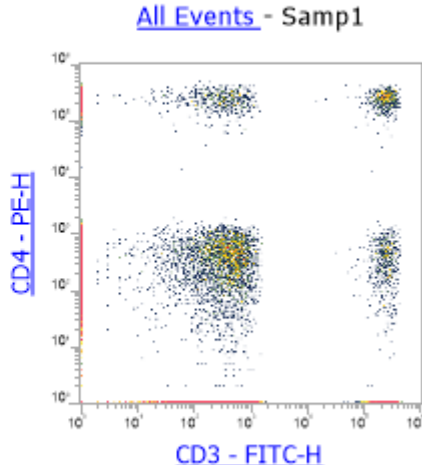
Tahoma A

**Plot Font Style**

**B**  
  *I*  
  U  
 10

**Axis Font Style**

**B**  
  *I*  
  U  
 9

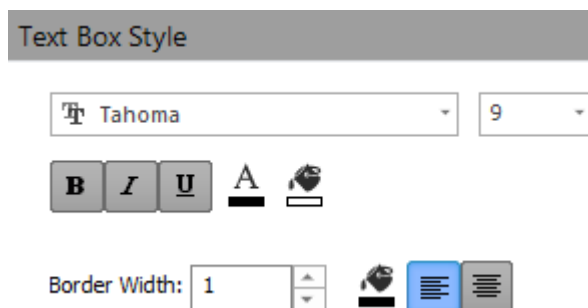


- Selecting the relevant **Include Name in Plot Title** options enables you to include the **Parent Gate**, **Experiment**, **Group**, and/or **Sample** name in the default plot title. The software appends the selections to the default plot title, separated by hyphens.
- Selecting the relevant **Include Name in Axis Label** options enables you to include the **Channel Name**, **Target** name, and/or **Label** name in the default axis labels. The software appends the selections to the default axis labels, separated by hyphens. By default, the **Target** and **Label** names are selected.
- **Font** options let you set the **Font** and **Font Color** for the plot title and axes labels. **Font** dropdown menu includes all fonts that are installed in the system. The default font is **Tahoma**. **Font Color** picker enables you to define the default font color of the plot title. The software uses the standard color picker dialog. The default color is **black**.
- **Plot Font Style** lets you select a font size from the dropdown menu, or enter a number from **6** to **72**. The default plot font size is **10 pt**.
- **Axis Font Style** lets select a font size from the dropdown menu, or directly enter a number. The default size in **9 pt**.
- You can select the **bold**, **italics**, and/or **underline** options to format the plot and axis fonts. Selecting one choice deselects the other.
- **Preview** section shows an example plot that reflects the current selections made in the **Fonts and Styles** tab. The preview is updated as the selections are changed.



## Text Box Style

**Text Box Style** section enables you to modify the default settings for text boxes.

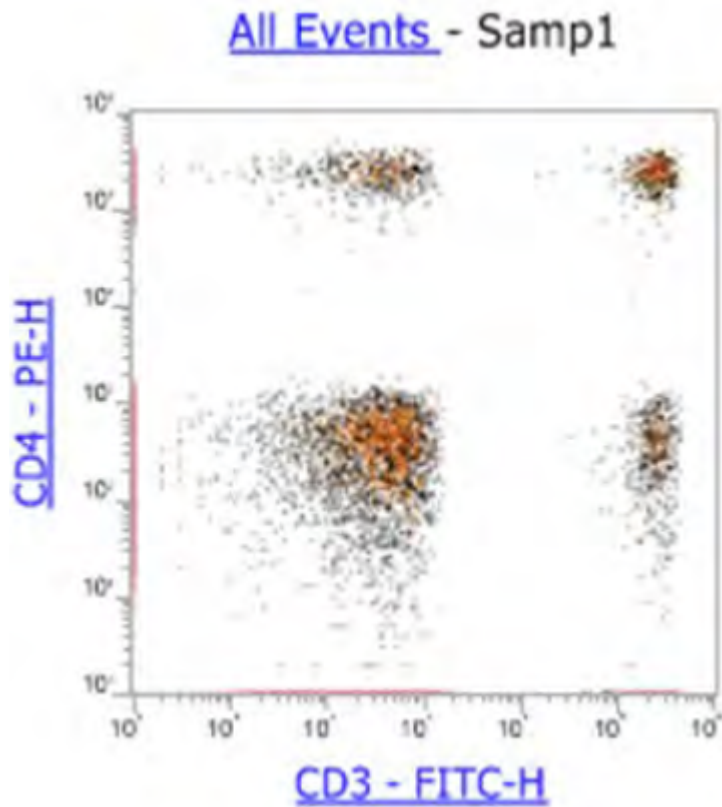


- You can select a font type from the **Font** dropdown menu. The dropdown menu includes all fonts that are installed in the system. The default font is **Tahoma**.
- In the **Font Size** field, you can select a font size from the dropdown menu, or directly enter a number. The default size is **9 pt**.
- You can select the **bold**, **italics**, and/or **underline** options to format the font. Selecting one choice deselects the other.
- Using the **Font Color** picker, you can define the default font color of the plot title. The software uses the standard color picker dialog. The default color is **black**.
- You can select to **left-align** or **center** the text. By default, text is **left-aligned**.
- You can select a **Border Width** for the text box, or enter an integer from **1** to **5**.
- You can select a **Border Color** and **Fill Color** using the appropriate color picker control. By default, border color is **black** and the text box is not filled.
- **Preview** section shows an example text box that reflects the current selections made in the **Fonts and Styles** tab. The preview is updated as the selections are changed.

## Previews

Previews section displays an example plot and an example text box reflecting the current selections made in the Fonts and Styles tab.

The previews are updated as the selections are changed.



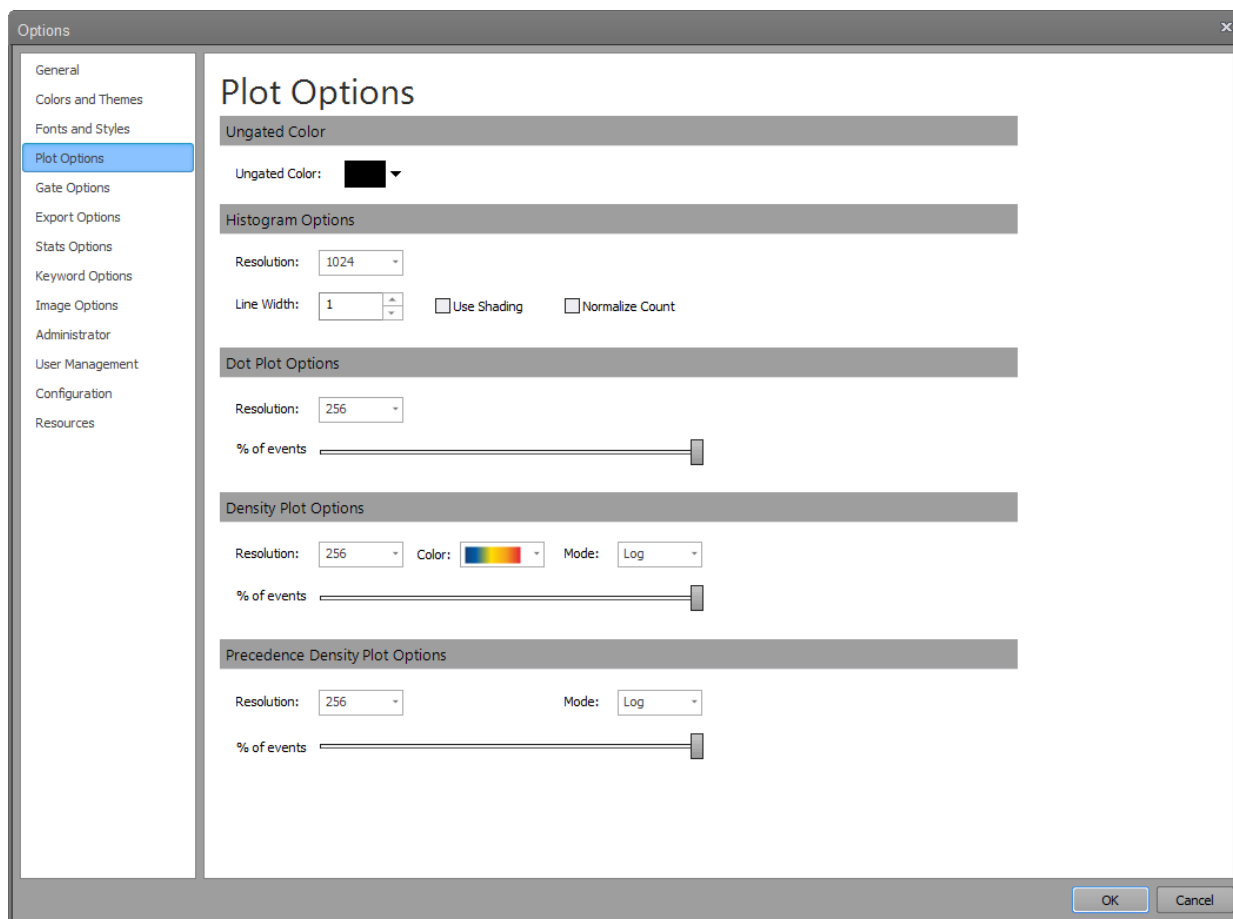
Text Box Style

# Plot Options

## Overview

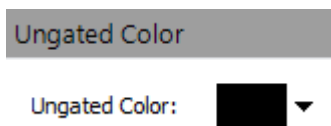
**Plot Options** is used to define display options for ungated plots (**Ungated Color**) and for each type of plot: **Histogram Options**, **Dot Plot Options**, **Density Plot Options**, and **Precedence Density Plot Options**.

Changes made in the Histogram plot, Dot plot, Density plot, and Precedence Density plot sections are only applied to newly created plots.



## Ungated Color

- By default, **Ungated Color** for all plot types is black.



- To modify the default color for ungated plots, click the **Ungated Color** dropdown button, which opens the standard color picker dialog.

- When a color is selected from the color picker dialog, the color displayed in the **Ungated Color** dropdown is updated and the color picker dialog is closed.
- Changes made to **Ungated Color** are applied immediately when **OK** is clicked and affect all past, present and future instances.

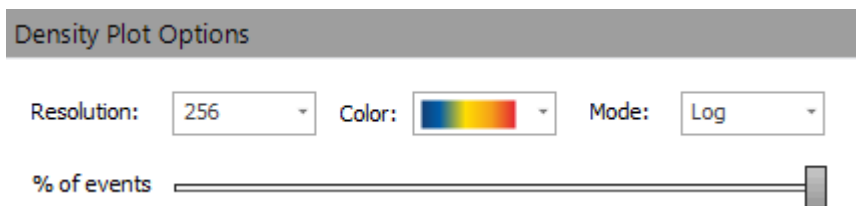
## Histogram Options

- **Resolution** enables you to modify the default resolution of Histogram plots. Available options are: **64, 128, 256, 512, and 1024**. The default resolution for Histograms is **1024**.
- **Line Width** enables you to modify the line thickness of Histogram plots. The minimum line thickness is **1 pixel** and the maximum is **5 pixels**. The default line thickness for Histograms is set at **1 pixel**.
- **Use Shading**, when selected, shades the area under the curve in the Histogram using the same color as the line color at 64% opacity. By default, **Use Shading** is deselected.
- **Normalize Histogram**, when selected, normalizes all new Histograms to the maximum peak. By default, **Normalize Histogram** is deselected.

## Dot Plot Options

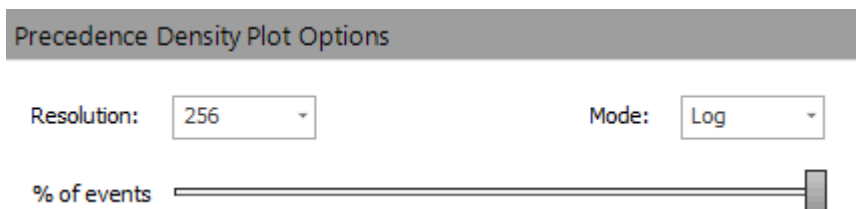
- **Resolution** enables you to modify the default resolution of Dot plots. Available options are: **64 × 64, 128 × 128, 256 × 256, 512 × 512, and 1024 × 1024**. The default resolution for Dot plots is **256 × 256**.
- **Density(%)** slider modifies the percentage of events displayed in a Dot plot from **0% to 100%** (left to right). The sample data are taken from the entire data file. By default, Dot plots show **100%** of events.

## Density Plot Options



- **Resolution** enables you to modify the default resolution of Density plots. Available options are: **64 × 64**, **128 × 128**, **256 × 256**, **512 × 512**, and **1024 × 1024**. The default resolution for Density plots is **256 × 256**.
- **Color** enables you to modify the default color setting for Density plots, which do not use the **Ungated Color** setting. The default color scheme for a Density plot is set to the first gradient available in the **Color** dropdown menu.
- **Mode** lets you change the **binning mode** for Density plots. Available options are **Log** and **Linear** modes. By default, the binning mode is set to **Log**.
- **Density(%)** slider modifies the percentage of events displayed in a Density plot from **0%** to **100%** (left to right). The sample data are taken from the entire data file. By default, Density plots show **100%** of events.

## Precedence Density Plot Options



- **Resolution** enables you to modify the default resolution of Precedence Density plots. Available options are: **64 × 64**, **128 × 128**, **256 × 256**, **512 × 512**, and **1024 × 1024**. The default resolution for Precedence Density plots is **256 × 256**.
- **Mode** lets you change the **binning mode** for Precedence Density plots. Available options are **Log** and **Linear** modes. By default, the binning mode is set to **Log**.
- **Density(%)** slider modifies the percentage of events displayed in a Precedence Density plot from **0%** to **100%** (left to right). The sample data are taken from the entire data file. By default, Precedence Density plots show **100%** of events.

---

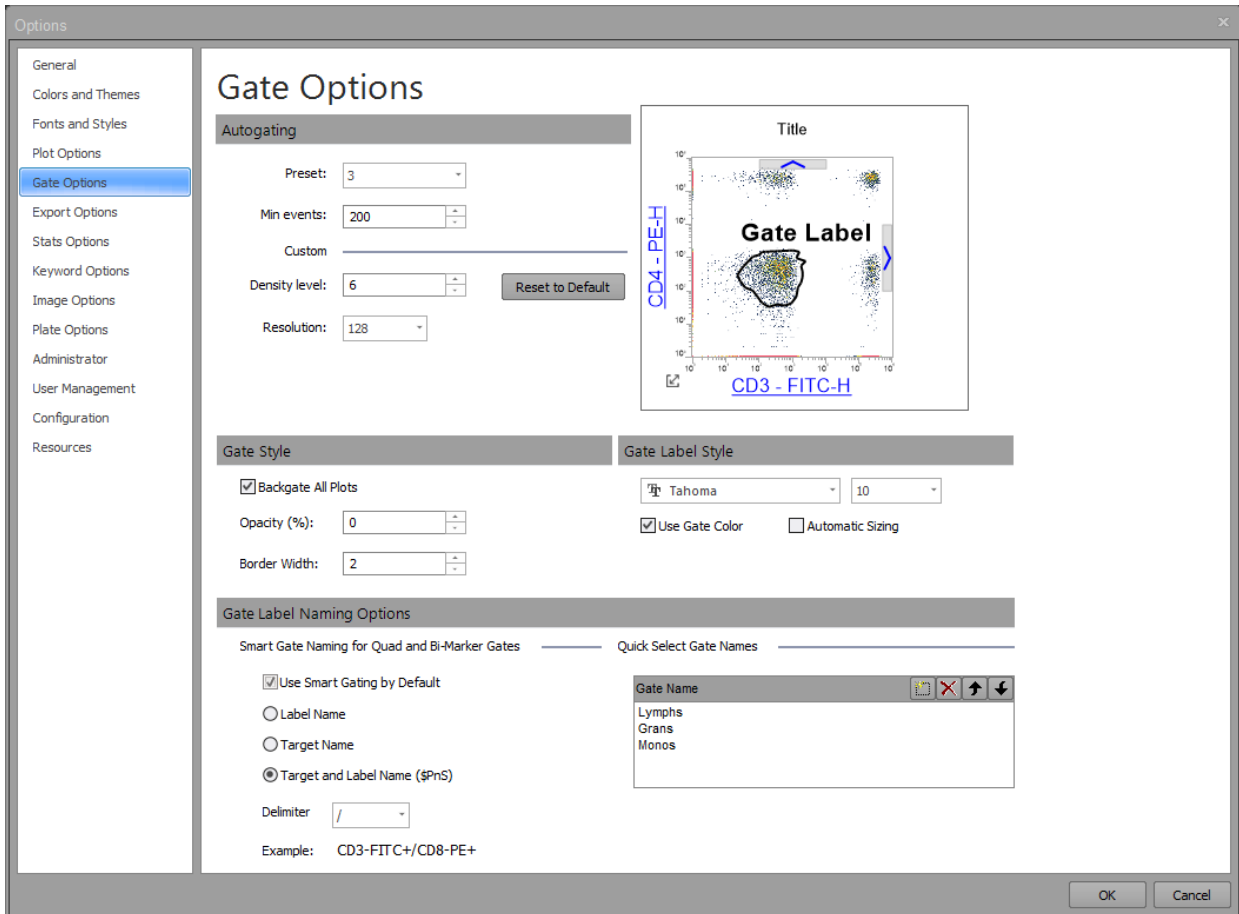
**Note:** The **Linear** mode bins data by assigning a color index for each density pixel linearly such that each increase is determined by dividing the range ( $Z_{\max} - Z_{\min}$ ) by the number of color steps. The **Log** mode bins data by assigning a color index for each density pixel logarithmically such that each increase is determined by dividing the logarithmic range ( $\log Z_{\max} - \log Z_{\min}$ ) by the number of color steps. The index is then determined as  $\text{int}((\log Z_{\text{val}} - \log Z_{\min}) / \text{increase})$ .

---

# Gate Options

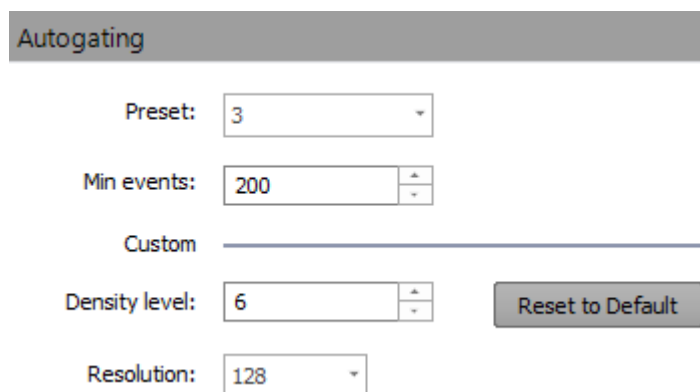
## Overview

**Gate Options** enables you to define the default settings for **Autogating**, **Gate Style**, and **Gate Label Style**, and for **Smart Gate Naming** and **Quick Select Gate Names** options.



## Autogating

**Autogating** enables you to set default autogating preferences, which include default preset, the number of events that must be contained in a gate, and custom settings for setting the density level and plot resolution.



The **Autogating** section contains the **Preset**, **Min events**, **Density level**, and **Resolution** controls, and the **Reset to Default** button (“Preset” on page 657).

---

**Note:** The autogating preferences that are defined in **Gate Options dialog** are only applied when creating new autogates. Changes to these settings do not affect autogate settings of existing gates. To change the settings of existing autogates, use the **Customize panel ▶ Gate options** (see “Customize gate options” on page 447).

---

### Preset

**Preset** dropdown enables you to set the default **Preset** for new autogates. A **Preset** consists of **Density level** and **Resolution** settings.

- The **Preset** dropdown has the following options:
  - **1** (Density level: 30, Resolution: 128)
  - **2** (Density level: 10, Resolution: 128)
  - **3** (Density level: 6, Resolution: 128)
  - **4** (Density level: 1, Resolution: 128)
  - **Custom** (By default, Density level: 5 and Resolution: 64, but you can change the custom settings using the **Density level** and **Resolution** controls)
- When a **Preset** is selected, the **Density level** and **Resolution** controls show the values that correspond to the **Preset** choice.
- By default, **Preset** is set at **3**.

## Min events

**Min events** sets the default minimum number of events that must be contained within a gate for the population to be autogated. This setting applies only to new autogates.

- If the **Min events** threshold is not reached and the gate is **ON**, the autogate remains in the **Autogate ON – [NO FIT]** mode.
- By default, **Min events** is set to **200**.
- The range of values for **Min events** is **1** to **9,999,999**.
- If you enter a **Min events** value that exceeds the upper numeric limit, then the value defaults to the maximum allowed number.

If the entered value for **Min events** is less than the lower limit, then the value defaults to the minimum allowed value.

## Density level

**Density level** sets the default **Density** value for the **Custom Preset** autogate option.

- **Density level** is populated with the value that corresponds to the selected **Preset** option (see Preset, above).
- The **Density** value for the **Custom Preset** listed above is the default value from a clean installation or for a new user account. The last used **Custom Preset** setting persists for the actively logged in user.
- The range for custom **Density** level is **1** to **100**.
- When the selected **Density level** is different from the selected **Preset**, the **Preset** dropdown reverts to **Custom**.

## Resolution

**Resolution** sets the default **Resolution** for the **Custom Preset** autogate option.

- **Resolution** field is populated with the value that corresponds to the selected **Preset** option (see Preset, above).
- The **Resolution** value for the **Custom Preset** listed above is the default value from a clean installation or for a new user account. The last used **Custom Preset** setting persists for the actively logged in user.
- The **Resolution** dropdown has the following options: **64, 128, 256, 512, 1024**.
- When the selected **Resolution** is different from the selected **Preset**, the **Preset** dropdown reverts to **Custom**.

## Reset to Default

**Reset to Default** sets the **Preset** to **3** and **Min events** to **200**.



## Gate Style

**Gate Style** options include settings for gate opacity, gate border width, and an option to add backgates on all plots.

Gate Style

Backgate All Plots

Opacity (%):

Border Width:

- **Opacity** enables you to set the gate opacity between **0%** and **75%**. By default, **Opacity** is set at **0%**.
- **Border Width** lets you set the gate border width between **1** and **5 pixels**. By default, **Border Width** is set to **2 pixels**.
- **Backgate All Plots** applies backgating to all **Oval**, **Rectangle**, and **Polygon gates** in all **Dot** and **Density plots**. Backgating paints the events contained within the gate on these plot types based on the color of the gate. By default, **Backgate All Plots** is selected.
- Clicking **OK** after changing the **Gate Style** options defines the new default settings for these options.
- Only newly created gates use the new default settings.

## Gate Label Style

**Gate Label Style** options include controls for gate label font and font size.

Gate Label Style

Tahoma

Use Gate Color  Automatic Sizing

- **Font** dropdown enables you to select a font type. The dropdown menu includes all fonts installed on the system. The default font is Tahoma.
- You can select a font size from the **Font Size** dropdown or enter a number from **6** to **72**.
- **Use Gate Color** checkbox sets the gate label color to match the gate color. By default, the checkbox is selected.  
When the checkbox deselected, the gate label color is black.
- **Automatic Sizing** checkbox enables the software to adjust the text size to fit in the plot. By default, the checkbox is not selected.
- Any changes made in this section are applied immediately when the dialog is closed and affect all past, present and future instances.

## Gate Label Naming

**Gate Label Naming Options** section includes controls for **Smart Gate Naming for Quad and Bi-Marker Gates** and **Quick Select Gate Names**.

Gate Label Naming Options

Smart Gate Naming for Quad and Bi-Marker Gates      Quick Select Gate Names

Use Smart Gating by Default

Label Name

Target Name

Target and Label Name (\$PnS)

Delimiter:

Example: CD3-FITC+/CD8-PE+

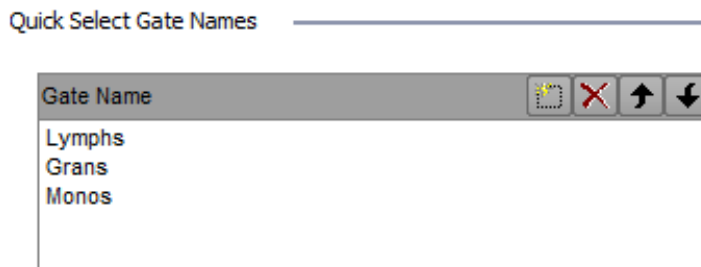
Gate Name

Lymphs  
Grans  
Monos

### Gate Label Naming Options

- **Use Smart Gate Naming** enables the software to automatically generate names for Quadrant gates. By default, the checkbox is selected.
- You can select the format for smart gate naming by clicking the appropriate control button. Available options are **Target Name**, **Label Name**, and **Target and Label Name (\$PnS)**. By default, smart gate naming format is set to **Target and Label name (\$PnS)**.
- For Quadrant gates, four gate names are generated (one for each quadrant).
- The **Delimiter** dropdown lets you to select a character to separate the names created for each quadrant of a Quadrant gate.
- An example of the concatenation of names and the delimiter is shown based on the selections made.
- Any changes made in this section are only applied to newly created gates.

## Quick Select Gate Names



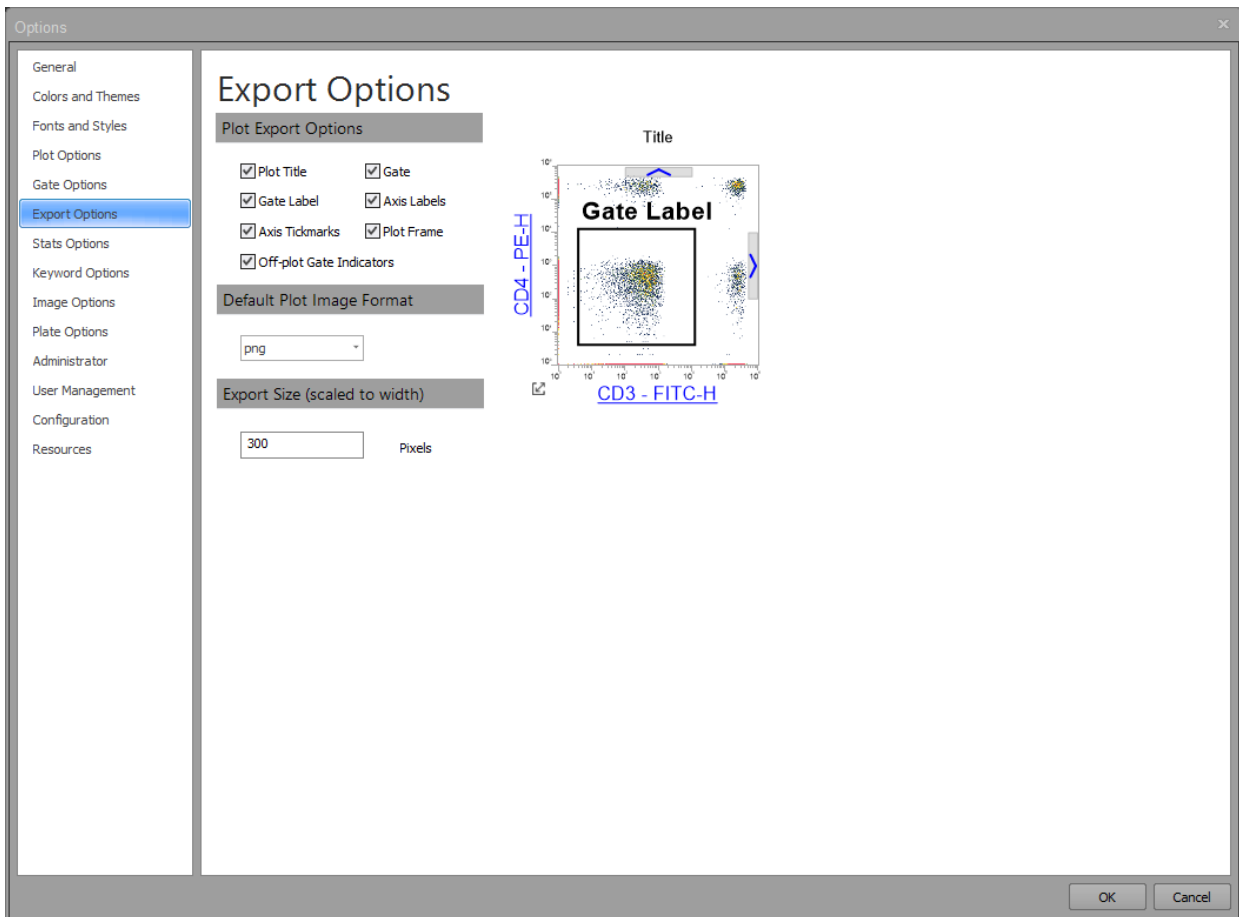
- **Quick Select Gate Names** lets you define the gate names available in the **Name** dropdown menu on the **Customize Gate** panel (“Name options” on page 459)
- By default, the **Quick Select Gate Names table** contains the **Lymphs**, **Grans**, and **Monos** options.
- You can add more names to the table by double-clicking on the whitespace of the list view or by clicking the **New** button in the table header.
- You can edit the names in the table by double-clicking on the name you want to edit.
- Gate names can be up to 50 characters long and can include any character except the following: \ / : \* < > | ?.
- Duplicate entries are not allowed.
- On validation, the software removes leading and trailing spaces, and converts consecutive spaces to single spaces.
- If you attempt to enter invalid characters, a warning dialog indicates the error condition, and the invalid characters do not appear in the text field.
- You can modify the sort order of the names in the table by selecting a row and using the **Up** and **Down** arrows in the table header.
- You can delete names from the table by selecting a row and using the **Delete** button in the table header.
- Changes made to entries in the table are available immediately. Gate names are not affected on established Workspaces.

# Export Options

## Overview

**Export Options** enables you to define the default plot content that is copied to the Windows™ clipboard using the **Copy** command (when a plot is pasted outside the application) or when the plot is saved using the **Save As** command. In addition, **Export Options** lets you specify the default plot image format used when saving plots.

Changes made in the **Export Options** tab take effect immediately and applied to newly exported objects.



## Plot Export Options

**Plot Export Options** let you select the default plot contents that are included when a plot is copied or saved. It consists of seven checkbox controls (**Plot Title**, **Gate**, **Gate Label**, **Axis Labels**, **Axis Tickmarks**, **Plot Frame**, **Off-plot Gate Indicators**), which are selected by default.

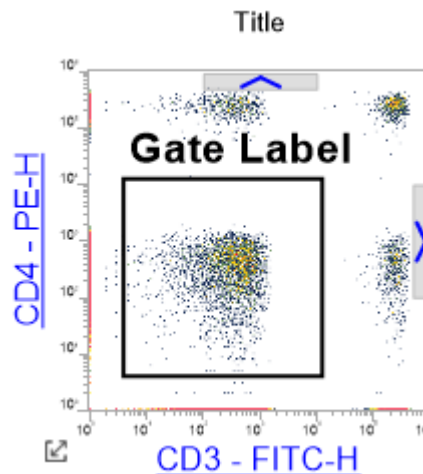
Plot Export Options

<input checked="" type="checkbox"/> Plot Title	<input checked="" type="checkbox"/> Gate
<input checked="" type="checkbox"/> Gate Label	<input checked="" type="checkbox"/> Axis Labels
<input checked="" type="checkbox"/> Axis Tickmarks	<input checked="" type="checkbox"/> Plot Frame
<input checked="" type="checkbox"/> Off-plot Gate Indicators	

- When the **Plot Export Options** checkboxes are selected or deselected, a **plot preview** shows what plot content is included when copying or saving a plot.

## Plot export preview

When the **Plot Export Options** checkboxes are selected, the **Plot export preview** updates to show which plot content will be included when copying or saving a plot.



## Default Plot Image Format

**Default Plot Image Format** option sets the default format used when saving a plot.

### Default Plot Image Format

png

- The **Plot Image Format** dropdown contains a list of available image formats. Available options are **PNG**, **GIF**, **JPG**, **TIF**, **BMP**, and **EMF**.
- By default, plot image format is set to **PNG**.

## Export Size (scaled to width)

**Export Size (scaled to width)** option enables you to set the **x-axis size in pixels** when exporting or copying objects.

### Export Size (scaled to width)

300      Pixels

- The x-axis pixel range for the export size is 100–5000 pixels. The default size is 400 pixels.
- When a size value greater than 5000 pixels is entered, the value defaults to 5000 pixels. When a value less than 100 pixels is entered, the value defaults to 100 pixels.

# Stats Options

## Overview

**Stats Options** let you define default display preferences for statistics boxes in the **Workspace** and **Results views**. The tab is divided into four sections: **Header**, **Statistics**, **Stats Box Style**, and **Decimal Settings**.

The screenshot shows the 'Stats Options' dialog box with the following settings:

- Header:**  Select All,  Plot Title,  Experiment,  Sample,  Time Recorded,  Group
- Statistics:**  Select All,  Plate,  Plot Title,  Volume,  X Median,  X %CV,  Experiment,  Gate,  Concentration,  Y Median,  Y %CY,  Group,  Autogate Status,  Percent Gated,  X Peak,  X rSD,  Sample,  X Parameter,  Percent Total,  Y Peak,  Y rSD,  Workspace,  Y Parameter,  X Mean,  X SD,  X r%CV,  Compensation Source,  Event Count,  Y Mean,  Y SD,  Y r%CV
- Concentration Units:** events/μL
- Stats Box Style:** Font: Tahoma, Size: 9, Bold: , Show Border:
- Decimal Settings:** Statistics Decimal Places: Concentration, 3

- Changes made in **Stats Options** are applied only to newly created statistics tables in the **Workspace** and newly created Experiments in the **Results view**, except **Decimal Settings**.
- Changes made to **Decimal Settings** are applied immediately when **OK** is clicked to close the dialog, and affect all past, present, and future instances.

## Header

**Header** options define the default format for the statistics box headers.

Header

Select All

Plot Title       Experiment       Group

Sample       Time Recorded

- You can select to show **Plot Title**, **Experiment**, **Group**, **Sample**, and **Time Recorded** in the header area of statistics boxes on the **Workspace**. By default, all options are selected.
- Selecting or deselecting **Select All** selects or deselects all other header checkboxes. Deselecting any of the checkboxes also deselects the **Select All** checkbox.

## Statistics

**Statistics** determine which default statistics are shown in statistics boxes on the **Workspace**.

Statistics

Select All

Plate       Plot Title       Volume       X Median       X %CV

Experiment       Gate       Concentration       Y Median       Y %CV

Group       Autogate Status       Percent Gated       X Peak       X rSD

Sample       X Parameter       Percent Total       Y Peak       Y rSD

Workspace       Y Parameter       X Mean       X SD       X r%CV

Compensation Source       Event Count       Y Mean       Y SD       Y r%CV

Concentration Units:

- You can select the default options you want to show in the statistics boxes by selecting the relevant **Statistics** checkbox. The following options are available:  
**Plate**, **Experiment**, **Group**, **Sample**, **Workspace**, **Compensation Source**, **Plot Title**, **Gate**, **X Parameter**, **Y Parameter**, **Event Count**, **Volume**, **Concentration**, **% Total**, **% Gated**, **X Mean**, **Y Mean**, **X Median**, **Y Median**, **X Mode**, **Y Mode**, **X SD**, **Y SD**, **X %CV**, **Y %CV**, **X rSD**, **Y rSD**, **X r%CV**, **Y r%CV**, and **Autogate Status**.  
You can change the statistics included in the statistics boxes at any time by selecting the statistics in the **Statistics Ribbon bar** when the **Statistics box** is selected.  
For more information about each option, see “Statistics tab” on page 103.
- By default, **Gate**, **X Parameter**, **Y Parameter**, **Event Count**, **% Gated**, **% Total**, and **Autogate Status** are selected.
- Default settings are applied to Experiment-level and Plot-level statistics boxes.
- Selecting or deselecting the **Select All** checkbox selects or deselects all statistics checkboxes. If any of the statistics checkboxes is deselected, the **Select All** checkbox is also deselected.



## Stats Box Style

**Stats Box Style** contains the formatting tools for the display of statistics boxes.

- You can select a **font type** from the **Font** dropdown menu. The **Font** dropdown menu includes all fonts installed on the system. The default font is **Tahoma**.
- In the **Font size** field, you can select a **font size** from the dropdown menu, or directly enter a number. The default size is **9 pt**.
- Using the **Font color** picker, you can define the default **font color** of the plot title. The software uses the standard color picker dialog. The default color is **black**.

## Decimal Settings

**Decimal Settings** define the default decimal places displayed for each statistic. It contains a combo box control with a list of statistics and a control to select the number of decimal places for each statistic.

- Clicking the **Statistics** combo box opens the **Statistics dropdown** list, which contains the following options:  
**All, Concentration, % Total, % Gated, Mean, Median, SD, %CV, rSD, r%CV, and Volume.**  
**Event Count** cannot be adjusted as it does not show any decimal places.
- When a statistic is selected from the list, the **Decimal Places** control updates to the default value assigned to the selected statistic.

The default number of decimals for each statistic is shown in the following table:

Statistic	Decimals
Concentration	3
% Total	3
% Gated	3
Mean	0
Median	0
SD	2
% CV	2
rSD	2

(continued)

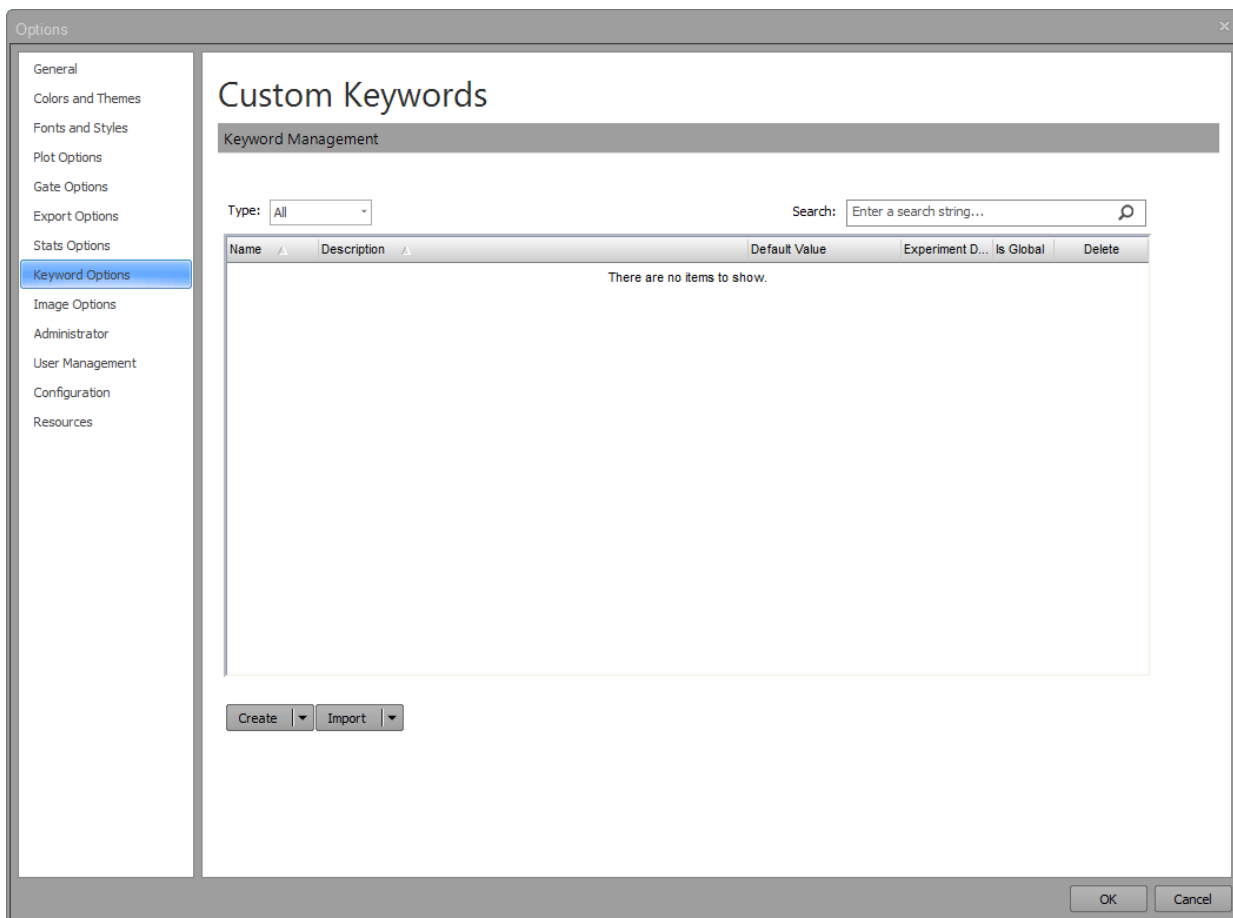
Statistic	Decimals
r%CV	2
Volume	0

- You can modify the number of decimal places displayed using the **Decimal Places** control. You can directly enter a number from **0** to **7**, or use the **up** and **down** buttons to modify the decimal places.  
Alternatively, you can use the **up** and **down** arrows on the keyboard, or the **mouse scroll wheel**.
- The text field only allows integer values. Entering a value greater than **7** defaults the value to **7**.
- The decimal places assigned to each statistic are used throughout the software, including the Workspace statistics tables, the Results tab values, the Report statistics tables, on-plot statistics for gates, and on the Heat Map legend.
- When **All** is selected from the **Statistic dropdown** list, the value input in the control applies to all statistics in the list.
- If the decimal places are different for each statistic, the control is left blank when **All** is selected.
- The number of decimal places specified here applies to all past, present, and future Experiments.
- The **Count** and **Mode** statistics only show integers without a decimal place.

# Keyword Options

## Overview

**Keyword Options** enable you to define custom keywords to provide more information about the Sample.



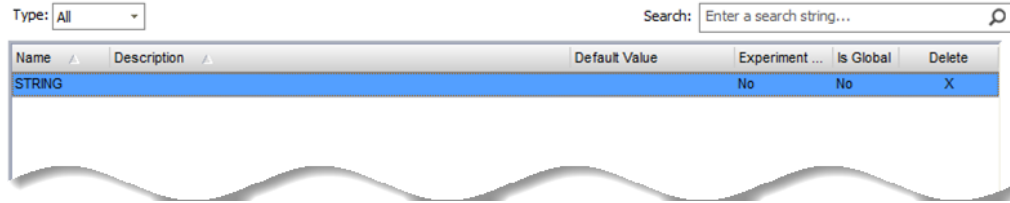
## Keyword Management

The **Keyword Options** tab contains a table displaying a list of all custom (i.e., user-defined) keywords and controls to create, edit, import and export keywords.

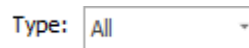
All users can create, edit and delete their own custom keywords and add them to their user list in **Keyword Options**.

## Keywords list

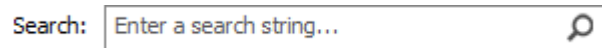
You can filter the keywords or search for existing keywords in the **keywords list**.



- Use the **Type** dropdown to filter existing keywords by type.

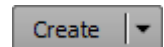


- Enter a search string into the **Search** field, then press **Enter** to search for existing keywords in the list.



## Create/Edit

**Create/Edit** split button enables you to create custom keywords and edit existing keywords with the **Create/Edit Keyword dialog** (“Create/Edit keyword dialog” on page 767).



- Click **Create** to open the **Create Keyword dialog**, then enter the name and description of the custom keyword you want to create.

- Select an existing keyword, then select **Edit** from the **Create/Edit dropdown** to open the **Edit Keyword dialog**.  
The **Edit Keyword dialog** is identical to the **Create Keyword dialog**, except the existing keyword values are populated.
- Only users with **Administrator** accounts can edit or delete a global keyword.  
Global keywords are displayed to all users, but the **Edit Keyword** option is disabled for non-administrator users.
- Edits made to a global keyword apply to the keyword master lists of all users, but they do **not** affect keywords that are associated with existing Experiments.
- To change the keyword included in an Experiment, use the **Edit** option for the keyword from the **Experiment Keyword dialog** (“Experiment Keywords dialog” on page 770) in the selected Experiment.

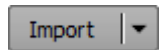
---

**Note:** After an Experiment is created, it gets its own set of keywords that persist with the Experiment. These keywords also persist through duplication, export/import, and as part of a Template.

---

## Export/Import

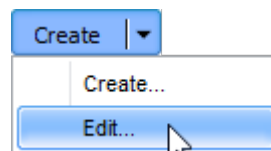
**Export/Import** split button enables you to export the selected keywords to a file, or to import keywords that were previously exported.



- Select an existing keyword, then click **Export** to export it to a file.  
The default file name for exported keywords is **AttuneKeywords** and the default file extension is **\*.akw**.
- If only a single keyword is selected, a single keyword is exported. To export multiple keywords, select multiple keywords (or the entire list), then click **Export**.
- Click **Import** to open the **File Open (Import) dialog** (“File Open (Import) dialog” on page 721), go to the appropriate folder, and import the desired keyword file (**\*akw**).  
If the imported keyword is a duplicate of an existing keyword, the duplicate keyword is appended by a numeric character (e.g., (1)).

## Set As Global

- **Set As Global** option in the **Create/Edit Keyword dialog** enables users with Administrator accounts to set the custom keywords in the Keyword options list to global. Global keywords are included in the Keyword Options for all user accounts.
- To set a custom keyword as global, select the keyword from the Keywords list, select **Edit** from the **Create/Edit dropdown**, then select **Set As Global** in the **Edit Keyword dialog**.



- The **Set As Global** option is only visible to users with **Administrator** accounts.

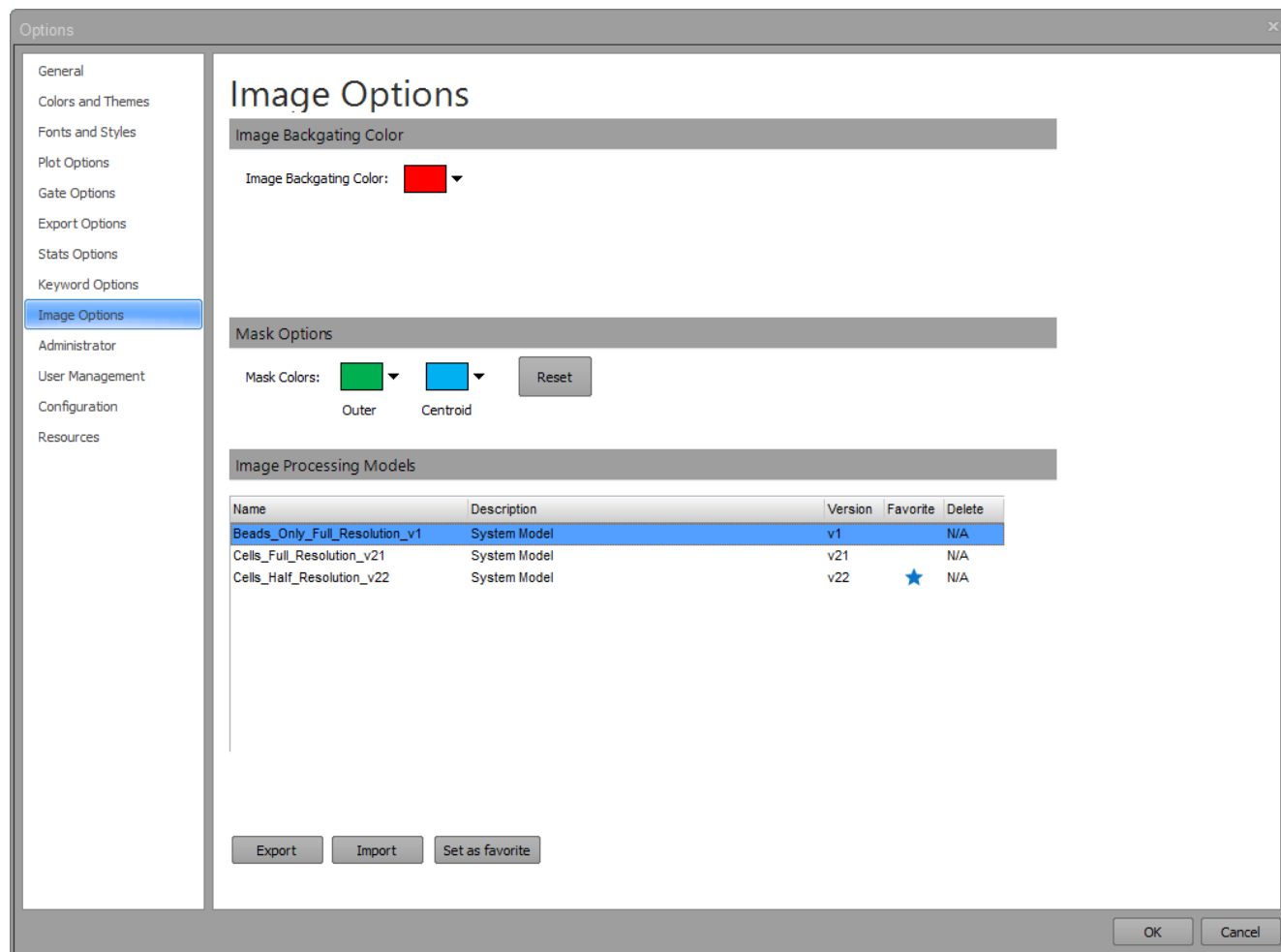
## Delete global keywords

- Only users with **Administrator** accounts can delete a global keyword.
- For non-administrator users, the **Keywords list** displays **N/A** in the **Delete** column for global keywords.
- Deleting a global keyword removes that keyword from the keyword master lists of all users, but it does not update existing Experiments.

## Image Options

**Image Options** tab contains the controls for **Image Backgating Color**, **Mask Options**, and **Image Processing Models**.

The **Image Options** tab is only available when using an Attune™ CytPix™ Flow Cytometer or when an Attune™ CytPix™ Experiment is active.



### Image Backgating Color

**Image Backgating Color** enables you to select the color with which the backgated imaged events are identified in the corresponding Workspace plots.

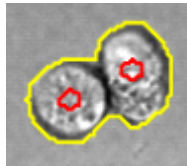
- You can modify the Image Backgating Color by clicking on the **color dropdown**, then selecting the new color from the standard **color picker dialog**.

When a color is selected from the **color picker dialog**, the color displayed in the **color dropdown** is updated and the **color picker dialog** is closed.

## Mask Options

**Mask Options** contain the controls to change the default **Outer** and **Centroid Mask Colors**.

- Image masks are binary representations of an image where an object is identified by the presence of a signal versus its absence. They provide a visual confirmation as to how the image processing “saw” the cell or the particle.
- The pixel coordinates of the mask contour are returned as outputs in the image processing results.
- The results include both the **outer masks** and the **centroid masks** (number of spots or cells detected in the Field of View). In the example below, the outer masks are depicted in yellow and the centroid masks in red.



- To change the default mask colors, select the desired **Mask Colors** for the **Outer** and **Centroid** masks, then click **OK** to close the **Options dialog**.
- To reset the default mask colors, click **Reset**, then click **OK** to close the **Options dialog**.

## Image Processing Models

**Image Processing Models** lists the available image processing models and lets you select the desired model as a **Favorite**. Image processing models are used in image processing operations. For a list and description of image processing models available with the Attune™ Cytometric Software, see Appendix B, “Attune™ Cytometric Software image processing parameters”.

---

**IMPORTANT!** Regardless of the size of the captured images in the experiment, images are processed at either full resolution (248 × 248 pixels) or at half resolution (124 × 124 pixels). Processing the images at half resolution decreases data footprint by 4-fold and improves the processing times by >4-fold. However, this comes with a potential trade off in accuracy.

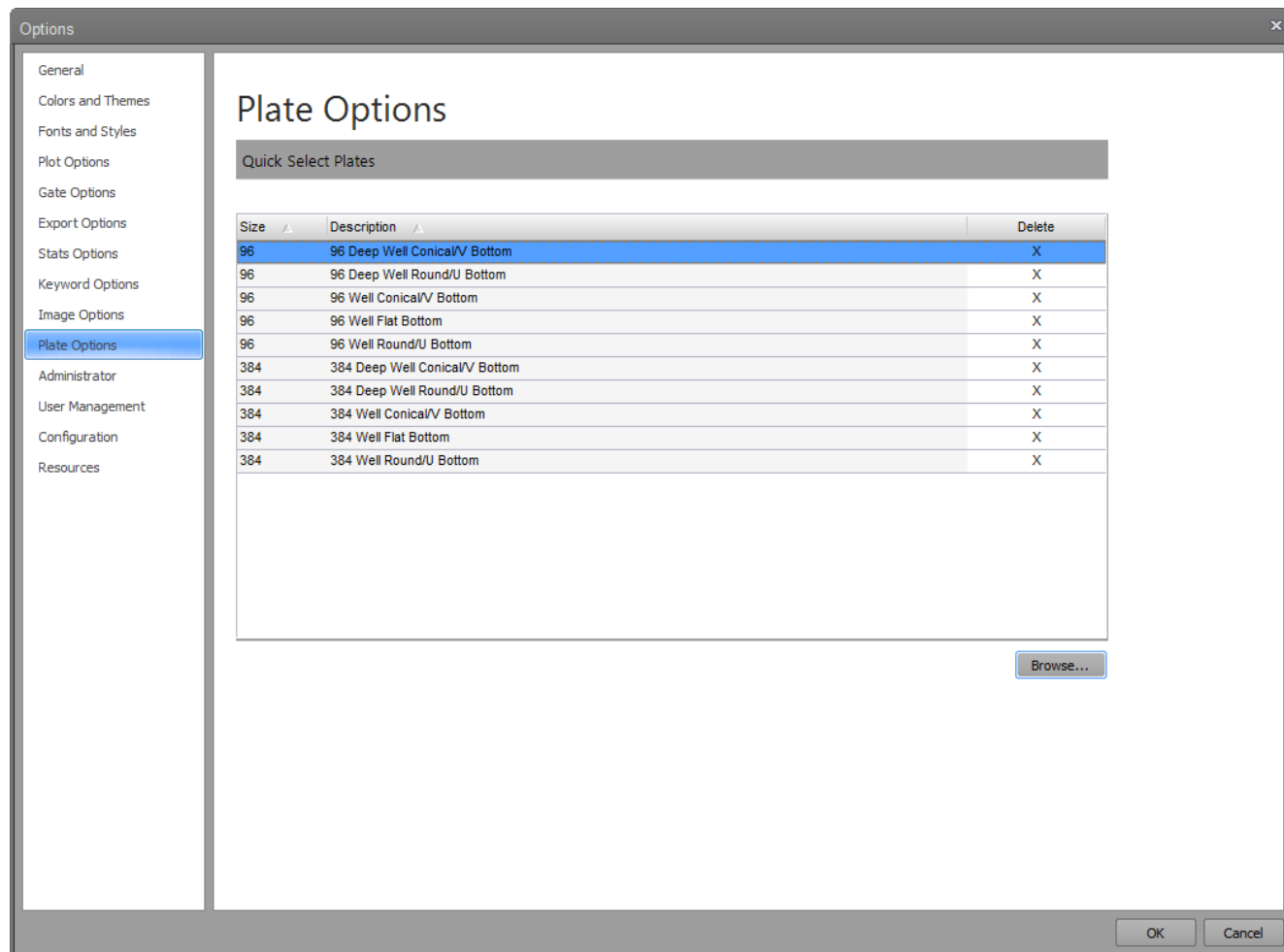
---

- To export an image processing model, select the desired model from the list, then click **Export** to open the **File Save (Export) dialog** (“File Save (Export) dialog” on page 715), go to the desired folder to save the image processing model file (with the file extension \*.dat), then click **Save**.
- To import an image processing model, click **Import** to open the **File Open (Import) dialog** (“File Open (Import) dialog” on page 721), go to the folder that contains the desired model, select the desired model (with the file extension \*.dat) from the list, then click **Open**.
- To set an image processing model as a **Favorite**, select the desired model from the list, then click **Set as favorite**.



## Plate Options

**Plate Options** tab contains the **Quick Select Plates** list, which enables you to designate selected plates as favorites. The selected plates become available in the **Plate Type** dropdown in the **New Experiment dialog** (“New Experiment dialog” on page 606) and in the **Customize Plate Experiment panel** (“Customize experiment (Plate or Tube) options” on page 466).



**Note:** The **Plate Options** dialog is available only when a CytKick™ Max™ Autosampler is connected to the Attune™ instrument.

- By default, the **Quick Select Plates** list includes the 10 defined plates available to all models. However, you can add extra plate definitions to the list using the **Select Plates** dialog (“Select Plates dialog” on page 774).  
To open the **Select Plates** dialog, click the **Browse** button.
- To sort the **Quick Select Plates** list, click the **column header**. The default sort order is plate size, followed by description.
- When the mouse moves over the **Delete** column, the cursor changes to the hand cursor, if the plate can be deleted. To remove a Plate from the list of favorites, click the **X** in the **Delete** column next to the plate you want to delete.

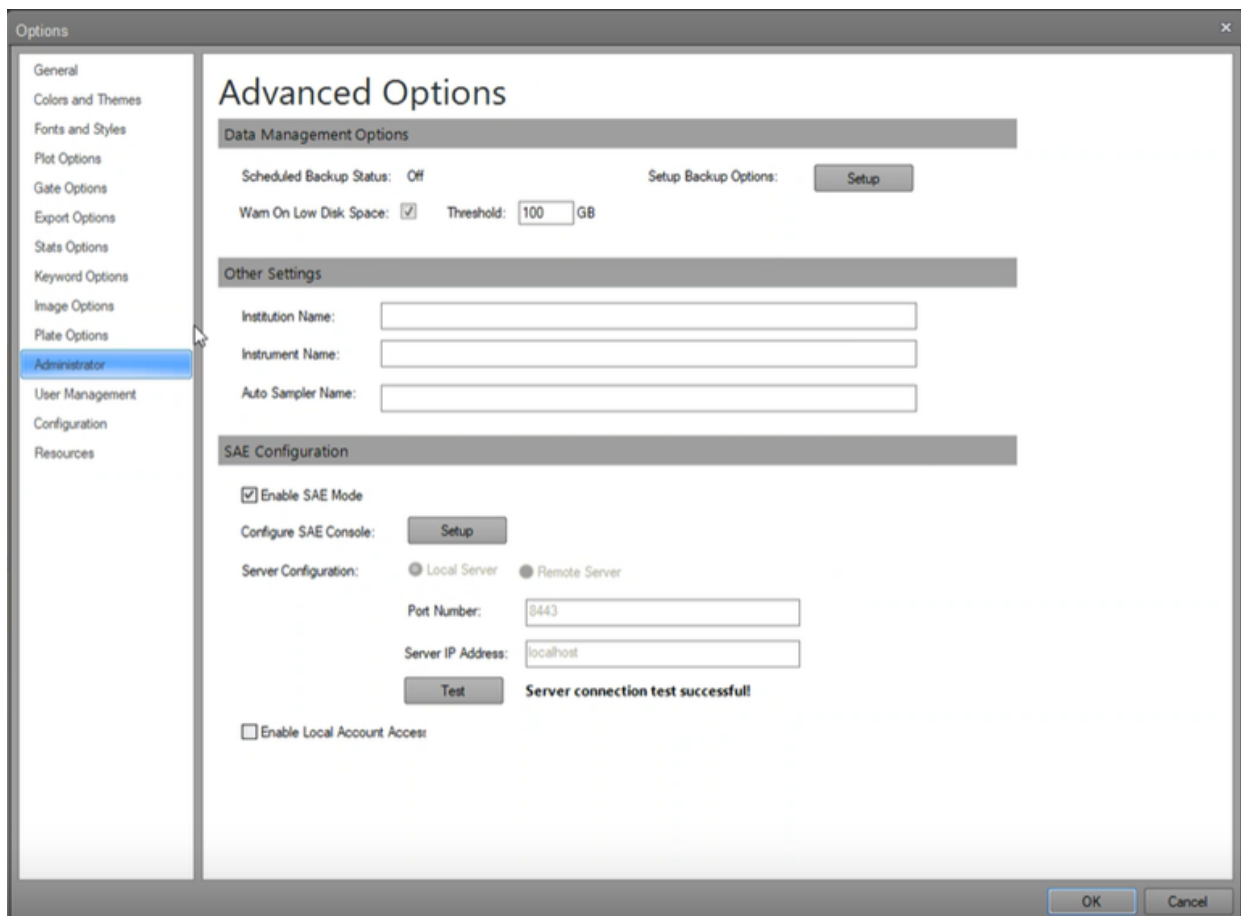
# Administrator

## Overview

The **Administrator** tab opens the **Advanced Options** controls, which enable users with Administrator account privileges to set:

- **Data Management Options:** Sets system level FCS backup policy (“Data Management Options” on page 677)
- **Other Settings:** Defines institution and instrument names (“Other Settings” on page 679)
- **SAE Configuration:** Enables the software to access the SAE (Security, Auditing, e–Signature) features (“SAE Configuration” on page 680)

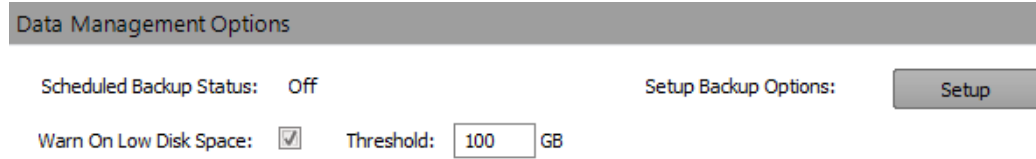
The **Administrator** tab is only visible to System Administrators, Administrators, and Service accounts.



**Note:** The **SAE Configuration Options** are only available if the SAE-specific DESkey device is present to enable access to the 21 CFR Part 11 features (see “SAE module” on page 37).

## Data Management Options

**Data Management Options** lets you to change backup and low disk space warning options, and to see the status of a scheduled backup.

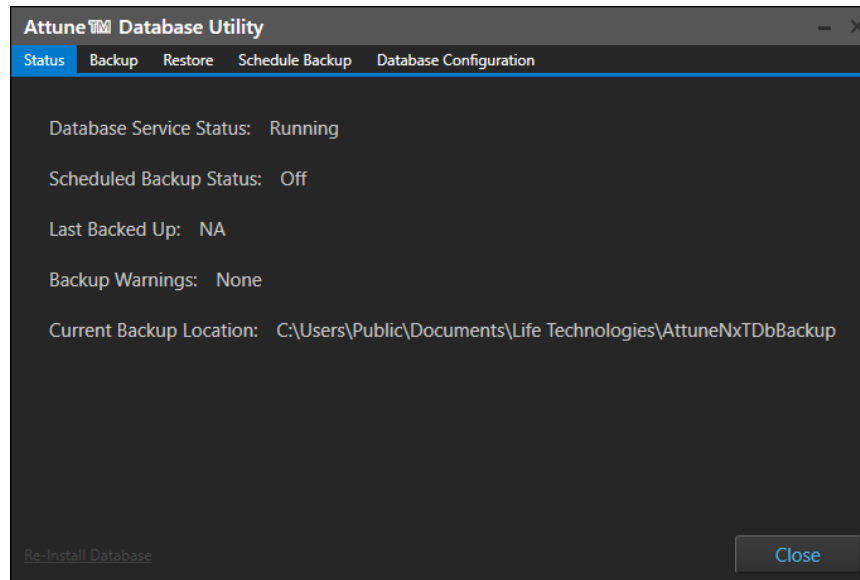


### Scheduled Backup Status

Shows whether the **scheduled backup** function is turned on or off.

### Setup Backup Options

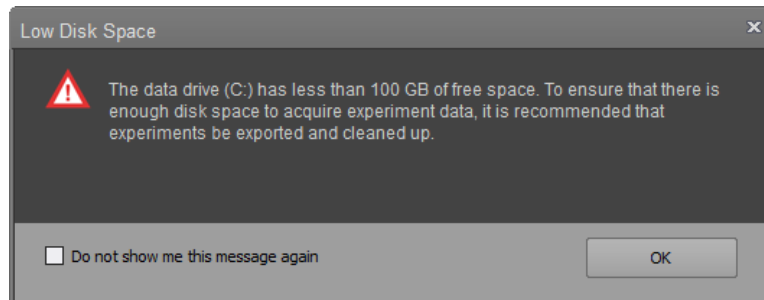
- Click **Setup** to open the Attune™ Database Utility (Chapter 29, “Attune™ Database Utility”).



- The database utility enables you to back up the data and database, to restore data and database, to schedule automated backups, or to reinstall the database.
- When the Attune™ CytPix™ Flow Cytometer is in use, the database utility enables you to select the Image Storage Path for captured images.
- By default, the captured images are stored at:  
C:\Users\Public\Public Documents\Life Technologies\AttuneCytPix\_Images

## Warn On Low Disk Space

Attune™ Cytometric Software displays a **Low Disk Space** warning when the available disk space on the primary disk drives falls below a preset **Threshold**.



- Users with Administrator or System Administrator accounts can change the low disk space warning options.
- To disable the **Low Disk Space** warning, deselect **Warn On Low Disk Space**.

Warn On Low Disk Space:  Threshold:  GB

- To change the threshold of available disk space on the primary disk drives below which the **Low Disk Space** warning is displayed, enter the desired value (in GB) in the **Threshold** field.

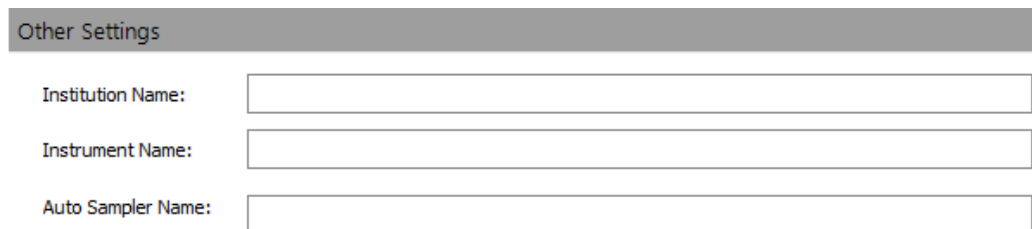
---

**Note:** By default, the **Threshold** is set to 100 GB. You can enter a threshold value of between 1 GB and 1000 GB.

---

## Other Settings

**Other Settings** includes settings for the **Institution Name**, **Instrument Name**, and **Auto Sampler Name**.



Other Settings

Institution Name:

Instrument Name:

Auto Sampler Name:

- You can enter up to 50 alpha-numeric characters in the **Institution Name**, **Instrument Name**, and **Auto Sampler Name** fields. The following characters are not allowed: \ / : \* < > |.
- If invalid characters are entered, a warning dialog indicates the error condition and the invalid characters will not appear in the text field.
- On validation, the software removes leading and trailing spaces.

### Institution Name

- To define the Institution Name for the system, type in the name into the **Institution Name** field.
- The Institution Name setting is used to populate the \$INST keyword in the FCS file TEXT segment (“Optional keywords” on page 829). The institution name is saved to each FCS file when entered in the **Institution Name** field.
- By default, the Institution Name is blank.

### Instrument Name

- To define the Instrument Name for the system, type in the name into the **Instrument Name** field.
- By default, the Instrument Name is blank.

### Auto Sampler Name

- To define the Auto Sampler Name for the system, type in the name into the **Auto Sampler Name** field.
- The Auto Sampler Name is blank by default.

## SAE Configuration

**SAE Configuration** options let you to set up the software to access the SAE (21 CFR Part 11) features.

- The **SAE Configuration** options are only available if the SAE-specific DESkey device is present to allow access to the 21 CFR Part 11 features (see “SAE module” on page 37).
- To configure the software for the **SAE mode**, the SAE server must be running.

### Enable SAE Mode

**Enable SAE Mode** option is used to enable and disable the **SAE mode** for the Attune™ Cytometric Software.

Enable SAE Mode

- To enable the **SAE mode**, check **Enable SAE Mode**.  
Only SAE user accounts with the Administrative role that have permission to **Configure Security and Auditing** can enable the SAE mode.  
For instructions about how to enable the **SAE mode**, see “Enable the SAE mode” on page 682.
- To disable the **SAE mode**, deselect **Enable SAE Mode**.  
Only local Attune™ users with Administrator or System Administrator account, or SAE user accounts with Administrator role can disable the **SAE mode**.  
For instructions about how to disable the **SAE mode** as an SAE user, see “Disable the SAE mode – SAE user” on page 684.  
For instructions about how to disable the **SAE mode** as a local Attune™ administrator or system administrator, see “Disable the SAE mode – local Attune™ administrator” on page 686.

---

**IMPORTANT!** When signing into the Attune™ Cytometric Software after the **SAE mode** is enabled, you need to use the SAE account username and password, and not the Attune™ Cytometric Software username and password.

Note that the data from local User and Advanced User accounts become inaccessible after the **SAE mode** is enabled, unless the **Enable Local Account Access** option is selected (see “Enable Local Account Access” on page 682). We recommend that you export all local User and Advanced User data before enabling the **SAE mode**.

---

## Configure SAE Administrator Console

**Configure SAE Administrator Console** lets you to configure SAE Console settings.

Configure SAE Console:

Setup

- To configure the SAE Administrator Console, click **Setup** to open the **SAE Administrator Console** window in the default web browser, then sign in to the **SAE Administrator Console**.
- You can create new SAE users, new SAE user roles, set security policies, and configure auditing and e-signature requirements in the **SAE Server Console** (see Appendix D, “SAE Administrator Console”).

## Server Configuration

**Server Configuration** enables you to configure the SAE server to be run on a local or remote server, and to set the Server IP Address and Port Number.

Server Configuration:

Local Server  Remote Server

Port Number:

8201

Server IP Address:

localhost

Test

- The **SAE Server Configuration** options are only enabled if the **SAE Mode** is disabled (i.e., **Enable SAE Mode** is unchecked; see “Disable the SAE mode – SAE user” on page 684).
- For both local and remote servers, the **Port Number** is set to 8201 by default and the **Server IP Address** is set to “localhost”.  
You can set the **Port Number** to 1025–49151.
- When **Local Server** option is selected, only the **Port Number** is editable. In this case, the **Server IP Address** is set to read-only and displays “localhost”.  
When the **Remote Server** option is selected, both the **Server IP Address** and **Port Number** can be changed.
- **Test** button enables you to test the connection to the **Remote Server**.  
When **Test** is clicked, the software attempts to connect to the **Remote Server** and performs a time check between the client and the server. For a successful connection, the time difference between the client and the server must be less than 5 minutes.
- If the test is successful, the “Server connection test successful!” message is displayed.

Server Configuration:

Local Server  Remote Server

Port Number:

8201

Server IP Address:

localhost

Test

**Server connection test successful!**

## Enable Local Account Access

**Enable Local Account Access**, when selected, gives permission to local User and Advanced User accounts to access their accounts when SAE mode is enabled.

Enable Local Account Access

- The local Administrator or SAE Administrator can select **Enable Local Account Access** at any point of usage.
- When **Enable Local Account Access** is selected, local accounts can access the Attune™ Cytometric Software, but **SAE Auditing** and **SAE e–Signature Control** will not occur for the local accounts.
- Disabling the **Enable Local Account Access** option does not prevent the local Administrator from accessing the Attune™ Cytometric Software.

## Rules for enabling the SAE mode

1. Only SAE User accounts with Administrative privilege that have permission to “Configure Security and Auditing” can enable the SAE mode.
2. Enabling the SAE mode requires a connection to the SAE server.
3. After the user provides credentials to enable the SAE mode, the following take place behind the scenes:
  - a. An authentication request for the administrator is sent to the SAE server.
  - b. If the provided credentials are valid and the specified SAE user account is authorized to enable the SAE mode, the server returns a success status code.
  - c. Upon successful administrator authentication, an SAE user session for the specified administrator account is initiated and all relevant security, audit, and e–Signature™ configurations take effect.

## Enable the SAE mode

1. To enable the **SAE mode**, check **Enable SAE Mode** (“SAE Configuration” on page 680).
2. When the **Enable SAE Mode** checkbox is checked, the software prompts you to perform a test with the proper IP address and port number.
3. Click **OK** to close the warning dialog, then enter the **Port Number** and **Server IP Address**.

Server Configuration:     Local Server     Remote Server

Port Number:   

Server IP Address:



- Click **Test** to test the connection to the server.

The software attempts to connect to the server and performs a time check between the client and the server. For a successful connection, the time difference between the client and the server must be less than 5 minutes.

If the test is successful, the “Server connection test successful!” message is displayed.

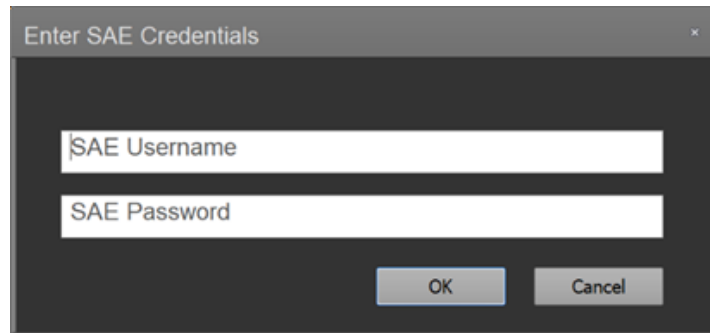
Server Configuration:       Local Server       Remote Server

Port Number:     

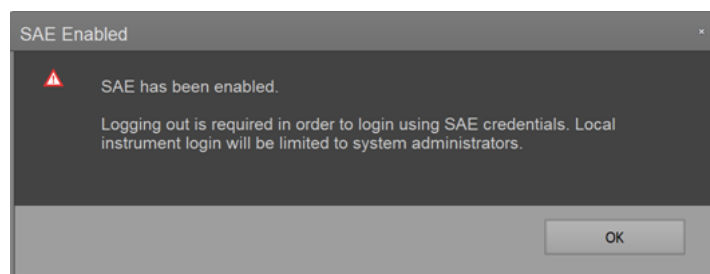
Server IP Address:     

     **Server connection test successful!**

- If the server can be reached and the test is successful, the **Enter SAE Credentials** dialog prompts you to enter the SAE credentials.



- If the credentials are correct, the **SAE Enabled** dialog is displayed.



If the server check fails, **Unable to connect to authentication server** dialog is displayed and the **Enable SAE Mode** is deselected.

- When the **SAE mode** is enabled, the Attune™ Cytometric Software displays the **SAE mode sign in** option (see “Sign in – SAE mode” on page 38).



All relevant security, audit, and e-signature configurations take effect immediately after enabling the SAE mode, and two action records appear in the SAE Administrator Console: “Enable Security” and “Login Success”.

### After the SAE mode is enabled

1. Only **SAE user** accounts can access the Attune™ instrument to perform runs (except **Administrator**, **System Administrator**, and **Service** users).
2. Users cannot sign in to the instrument via the local instrument profiles that were in use previously:
  - a. The local instrument profiles and their associated data are not deleted via this action of enabling SAE.
  - b. The action of disabling SAE enables these local instrument profiles to be accessible again.

---

**IMPORTANT!** When the application is configured to the **SAE mode**, you can sign in to the local Attune™ Cytometric Software account only if you have an **Administrator**, **System Administrator**, or **Service** account. **User** and **Advanced User** accounts are not allowed to sign in to the local account after SAE is enabled.

---

## Disable the SAE mode – SAE user

### Rules for disabling the SAE mode as an SAE user

1. Only SAE user accounts that have **SAE Administrator** privileges can disable the **SAE mode**.
2. Disabling the **SAE mode** requires a connection to the SAE server.

---

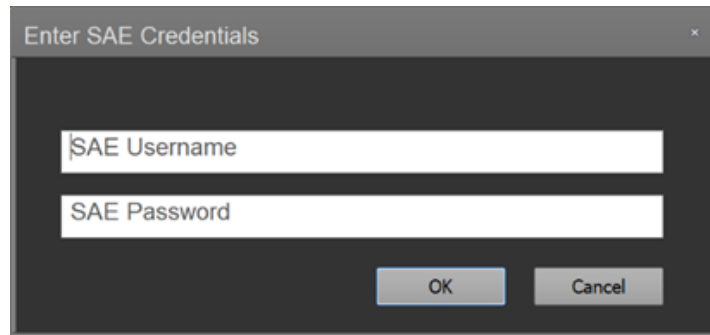
**Note:** Local Attune™ users with **Administrator** or **System Administrator** accounts can disable the **SAE mode** regardless of server status (“Disable the SAE mode – local Attune™ administrator” on page 686).

---

3. After the SAE user provides credentials to disable the **SAE mode**, the following take place behind the scenes:
  - a. An authentication request for the **SAE Administrator** role is sent to the SAE server.
  - b. If the provided credentials are valid and the specified SAE user account is authorized to disable the **SAE mode**, the server returns a success status code.
  - c. After successful administrator authentication, the instrument logs out of the current SAE account, and all relevant security, audit, and e-Signature configuration no longer affect the instrument.

## Disable the SAE mode

1. To disable the **SAE mode**, deselect **Enable SAE Mode** (“SAE Configuration” on page 680).
2. If the server is online and can be reached, the **SAE Credentials dialog** prompts you to enter the SAE credentials.



The image shows a dialog box titled "Enter SAE Credentials". It contains two text input fields, one labeled "SAE Username" and one labeled "SAE Password". At the bottom of the dialog, there are two buttons: "OK" and "Cancel".

3. The credentials are checked on the SAE server and only users with the permission to “Configure Security and Auditing” can disable **SAE mode**.  
If an error occurs when entering credentials, an error message is displayed in the **SAE Credentials dialog**.
4. If the entered credentials are correct, the **SAE Credentials dialog** closes and the SAE user is logged out of the application. The login screen displays the non-SAE mode sign in options (see “Sign in – Standard mode” “Sign in – standard mode” on page 38).



The image shows a login screen with two text input fields, one labeled "Username" and one labeled "Password". Below the fields, there is a link labeled "Forgot Password?" and a button labeled "Sign in".

All relevant security, audit, and e-Signature configurations are disabled immediately after disabling the **SAE mode**, and two action records appear in SAE Administrator Console: “Disable Security” and “Logout”.

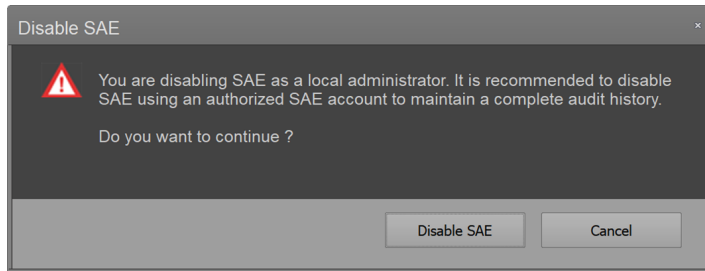
## Disable the SAE mode – local Attune™ administrator

### Rules for disabling the SAE mode as a local Attune™ user

1. Only local Attune™ users with **Administrator** or **System Administrator** accounts can disable the **SAE mode**.
2. Local Attune™ users with **Administrator** or **System Administrator** accounts can disable the **SAE mode** regardless of server status.

### Disable the SAE mode as a local Attune™ user

1. To disable the **SAE mode**, deselect **Enable SAE Mode**.
2. When **Enable SAE Mode** is deselected, a warning dialog is displayed to confirm that you want to disable the SAE.



3. Click **Disable SAE** to disable the account or click **Cancel** to close the dialog without disabling the **SAE mode**.
4. When SAE is disabled by a local Attune™ user, the “Disable Security” event is audited on the SAE Server in **Audit History ► Action Records** page.

---

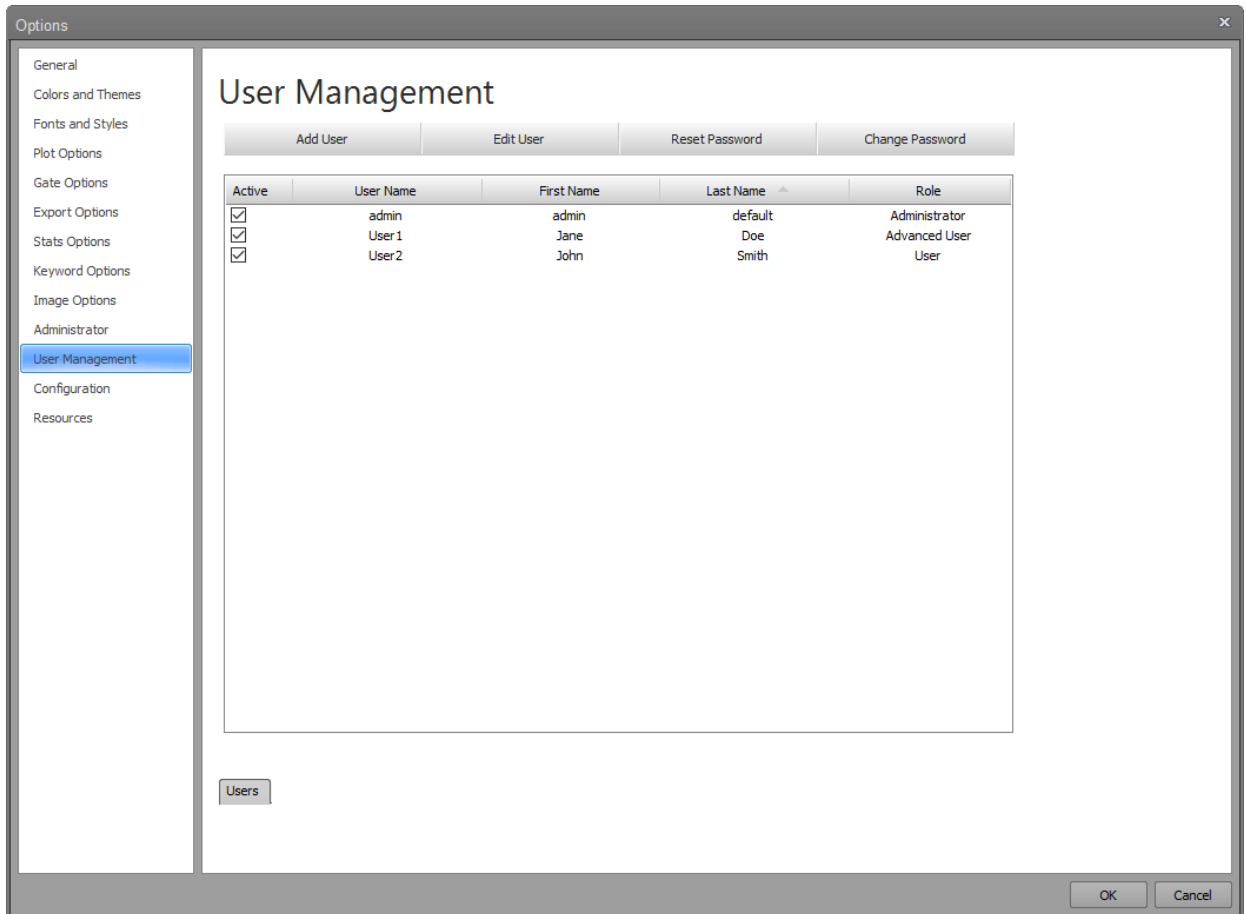
**Note:** To maintain a complete audit history, we recommend that you disable the **SAE mode** using an authorized SAE account (“Disable the SAE mode – SAE user” on page 684).

---

# User Management

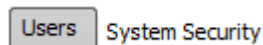
## Overview

**User Management** enables authorized users to add, edit, and view local Attune™ user accounts and manage user account passwords.



**Note:** When the **SAE mode** is enabled (see “Enable the SAE mode” on page 682), user management is controlled by the SAE Administrator Console (see Appendix D, “SAE Administrator Console”). For SAE specific permissions, see “SAE account permissions” on page 887.

- The tabs at the bottom of the **User Management** indicate the menu selected.



**Note:** For all account types except for the **System Administrator**, only the **Users** tab (“Users tab” on page 688) is visible. **System Administrators** have access to define the **System Security** (“System Security tab” on page 692) setting in the **System Security** tab.

- The **User Management** tab provides users with different account types the privileges described in “Default local Attune™ accounts” on page 42.

- For **Users**, the **Users** tab (“Users tab” on page 688) contains the controls for editing and viewing user’s accounts and for changing the user account passwords.  
For **Administrators**, the **Users** tab (“Users tab” on page 688) contains controls for adding users, editing and viewing accounts, and resetting or changing passwords.  
For **System Administrators**, the **Users** tab (“Users tab” on page 688) contains controls for adding and editing user accounts and for resetting or changing passwords. The System Administrator can also delete user accounts from the Users tab.
- User privileges for different account types are described in “Account permissions” on page 44.
- The settings defined in the **User Management** tab are global and they are stored in and retrieved from the database.

## Users tab

**Users** tab enables authorized users to add and edit users and to reset or change passwords. The table of User Accounts below these controls displays the database-derived information for all users.

Any changes made in the **Users** tab take effect immediately and are not reversed if **Cancel** is clicked in the **Options dialog**.

Add User      Edit User <b>Reset Password</b> Change Password				
Active	User Name	First Name	Last Name ▲	Role
<input checked="" type="checkbox"/>	admin	admin	default	Administrator
<input checked="" type="checkbox"/>	User1	Jane	Doe	Advanced User
<input checked="" type="checkbox"/>	User2	John	Smith	User

Users

Figure 135 Users tab – Administrator account view

Add User						Edit User						Reset Password						Change Password					
Active	User Name	First Name	Last Name	Role	Delete																		
<input checked="" type="checkbox"/>	service	service	default	Service	[X]																		
<input checked="" type="checkbox"/>	admin	admin	default	Administrator	[X]																		
<input checked="" type="checkbox"/>	sysadmin	sysadmin	default	System Administrator	[X]																		
<input checked="" type="checkbox"/>	User 1	Jane	Doe	Advanced User	[X]																		
<input checked="" type="checkbox"/>	User 2	John	Smith	User	[X]																		

Users System Security

Figure 136 Users tab – System Administrator account view

## Add User

- **Add User** enables an authorized user to create a new user account. This button is enabled only for **Administrator** and **System Administrator** accounts.
- Clicking **Add User** opens the **Add User dialog** (“Create a user profile” on page 706).
- If **SAE mode** is enabled (“Enable the SAE mode” on page 682), new SAE user accounts can only be created using the SAE Administrator Console (see **Users tab** in “Users tab” on page 882).
- The **Add User** button is disabled if an SAE user is signed in.

## Edit User

- **Edit User** enables an authorized user to modify a user account profile or the current user to modify their user profile.
- The **Edit User** button is enabled for all local Attune™ users.
- **Administrator** and **System Administrator** accounts can edit any user account as described in “Edit a user profile” on page 708. Any user can edit their own account.
- Clicking **Edit User** opens the **Edit User dialog** (“Edit a user profile” on page 708).
- SAE users cannot edit their user profiles using the Attune™ software and must use the SAE Administrator Console (see **Users tab** in “Users tab” on page 882)).
- The **Edit User** button is disabled when signed in as an SAE user.

## Reset Password

- **Reset Password** enables an authorized user to reset the password for any user account.
- The **Reset Password** button is enabled only for local Attune™ **Administrator** and **System Administrator** accounts.  
SAE users who have the SAE permission to “Configure Security and Auditing” cannot reset user passwords using the Attune™ software and must use the SAE Administrator Console (see **Users tab** in “Users tab” on page 882).
- Clicking **Reset Password** opens the **Reset Password dialog** (“Reset a password” on page 711).

## Change Password

- **Change Password** enables the current users to change their own passwords. This button is enabled for all users.
- Clicking **Change Password** opens the **Change Password dialog** (“Change a password” on page 712).
- **Change Password** function is available for both SAE and local Attune™ users.

---

**Note:** For local Attune™ users, the password policy applied to new and changing passwords is based on the system security policy for user passwords set in the **System Security** tab of **User Management** options (see “System Security tab” on page 692).

For SAE users, the password policy applied to new and changing passwords is set based on the SAE security policy (see “System tab” on page 897 in Appendix D, “SAE Administrator Console”).

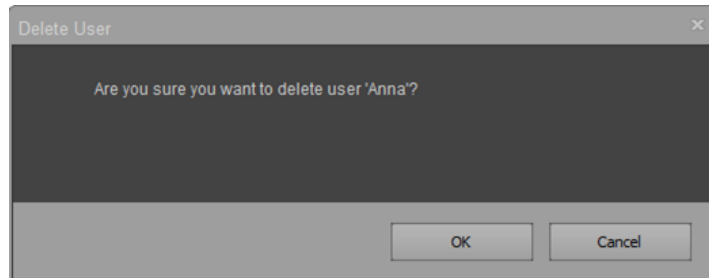
---

## User Accounts table

- The **User Accounts** table displays a list of all user accounts. It consists of columns for **Active** (i.e., **Account Status**), **User Name**, **First Name**, **Last Name**, **Role**.  
For **System Administrator** accounts, the table also has a **Delete** column with an **X** (i.e., **Delete**) button to remove specific user accounts.
- All selections in this table are single select. No multi selection is allowed.
- The table can be sorted by clicking on the column headers. When clicked, a sort indicator in the column heading specifies both the active sort column and the sort direction. Any single column can be used for sorting. Clicking on the same header reverses the sort direction.
- The default sort order is ascending by **Last Name**.
- Any user can see their own account listed in the User Accounts table. Only users with **Administrator** and **System Administrator** accounts can view other user accounts in the table.
- The **Active** (i.e., **Account Status**) column is only enabled for users with **Administrator** and **System Administrator** accounts to allow enabling and disabling of accounts. All other users can view the column, but the checkbox is disabled.
- When the **Active** checkbox is selected, the relevant account is enabled. When unchecked, the account is disabled.
- Disabling an account prevents a user from signing in. If a user with a disabled account attempts to sign in, the software displays a warning message to notify that the account has been disabled.
- The **Delete** option is only visible and enabled for the **System Administrator** account.



- When the **X** (i.e., **Delete**) button is clicked, the **Delete User dialog** prompts the user to confirm the deletion. Only a single user can be selected for deletion at a time.



- Clicking **OK** deletes the account, closes the **Delete User dialog**, and updates the **User Accounts** table.  
Clicking **Cancel** closes the **Delete User dialog** without deleting the account.
- When an account is deleted, it is removed from the list of user accounts visible to **Administrator** accounts.
- Only the **System Administrator** can view the full list of all users, where deleted user accounts are listed with the **Active** checkbox unselected and disabled, and the **Role** column set to **Deleted**.
- When an account is deleted, the account is also set to a **Deleted** status, and the **Active** checkbox is unselected and disabled. The account remains in the list.
- A deleted account relinquishes the username, allowing the username to be reused.
- SAE user accounts and local service accounts cannot be set to inactive or be deleted. If the **System Administrator** account attempts to delete or disable an SAE user account, a warning dialog is displayed.

## System Security tab

**System Security** tab lets the **System Administrator** define the system security policy for user passwords and password expiration criteria. This tab is only accessible to **System Administrator** accounts.

**User Names**

The length of the username field must be between 3 and 50 characters  
No spaces are allowed.

**User passwords**

Minimum password length:

User may not reuse the previous  passwords.

Password must contain a least one of the following characters:

Uppercase       Numeric  
 Lowercase       Non alpha-numeric (e.g !, # < \$, %)

**Password Expiration**

Password expires automatically after  days.  
Notify the user  days before expiration.

Lock out user after  failed attempts.  
Wait  minutes after locking out user to allow further login

Users **System Security**

## User passwords

**User passwords** section provides options to set the minimum password length, to limit the reuse of passwords, and to set password complexity requirements.

User passwords

Minimum password length:

User may not reuse the previous  passwords.

Password must contain a least one of the following characters:

Uppercase       Numeric

Lowercase       Non alpha-numeric (e.g !, # < \$, %)

- The default **Minimum password length** is 4 characters.
- The default for **User may not reuse the previous passwords** is 0, which lets users reuse their last password.
- The maximum allowed value for password reuse is 10, which prohibits the use of the previous 10 passwords. If a value greater than 10 is entered, the value defaults to 10.
- The **Password complexity requirements** are unchecked by default. When selected, passwords are required to have at least one character from the selected complexity options.
- **Uppercase** selection requires that at least one uppercase alphabetic character is used in a password. Uppercase characters are defined from European languages (A through Z, with diacritic marks, Greek and Cyrillic characters).
- **Lowercase** selection requires that at least one lowercase alphabetic character is used in a password. Lowercase characters are defined from European languages (a through z, sharp-s, with diacritic marks, Greek and Cyrillic characters).
- **Numeric** selection requires that at least one numeric character is used in a password. A numeric character is defined as a base 10 digit (0 through 9).
- **Non-alpha-numeric** selection requires that at least one non-alpha-numeric character is used in a password. The non-alpha-numeric characters are: ~ ! @ # \$ % ^ & \* \_ - + = ` | \ ( ) { } [ ] ; : " ' < > , . ? /.
- Any Unicode character that is categorized as an alphabetic character but is not uppercase or lowercase can also be used. This includes Unicode characters from Asian languages.

## Password expiration

**Password expiration** section is used to set a period to force users to change their passwords, to notify users before their password expires, and to lock out users after a defined number of failed login attempts.

**Password Expiration**

Password expires automatically after  days.

Notify the user  days before expiration.

Lock out user after  failed attempts.

Wait  minutes after locking out user to allow further login

- For all number fields, if you enter a number that is lower than the minimum allowed, the number is adjusted to the minimum when you elsewhere or when you press **Enter**.
- If you enter a number that is greater than the maximum allowed, the number is adjusted to the maximum when you elsewhere or when you press **Enter**.
- Only integers are allowed.
- The default setting for **password expiration** is 90 days and deselected.
- The minimum value for **password expiration** is 30 days and the maximum is 365 days.
- The default setting for **password reset notification** is 10 days before expiration and deselected.
- The minimum value for notification is 1 day and the maximum is 10 days before password expiration.
- The default setting for **password lockout** for failed login attempts is 5 failed attempts and selected.
- The minimum value for **password lockout** is 3 failed attempts and the maximum is 10 failed attempts.
- The default setting for **wait between failed login attempts** is 5 minutes and selected.
- The minimum value for **wait between failed login attempts** is 1 minute and the maximum is 15 minutes.

# Configuration

## Overview

*Configuration* options allow you to set the offline instrument configuration when working in analysis-only conditions, and to manage the bubble sensor and automation port settings.

Options

General  
Colors and Themes  
Fonts and Styles  
Plot Options  
Gate Options  
Export Options  
Stats Options  
Keyword Options  
Image Options  
Plate Options  
Administrator  
User Management  
Configuration  
Resources

### Access resources for Attune™ Cytometer Hardware

Hardware / Virtual Laser Configuration

Instrument Type: Attune CytPix Flow Cytometer

Attune Model: 0A48663 Attune CytPix Flow Cytometer (Lasers: BRVY)

Autosampler Model: CytKick™ Max Autosampler

Enable Bubble Sensor

Enabled

Automation Communication Settings

TCP/IP Communication Port: Valid ports are: 1025-49151 (Default is 8225)

Thermo Fisher Cloud Device Registration

Device Name:

Register Device

OK Cancel

Configuration options tab contains the following controls:

- *Hardware/Virtual Laser Configuration* (“Hardware/Virtual laser configuration” on page 696)
- *Enable Bubble Sensor* (“Enable bubble sensor” on page 697)
- *Automation Communication Settings* (“Automation communication settings” on page 697)
- *Connect Device Registration* (“Connect device registration” on page 698)

## Hardware/Virtual laser configuration

*Hardware/Virtual Laser Configuration* allows you select offline instrument type and model. This selection aids in the creation of Experiments when not connected to an Attune™ instrument by maintaining correct instrument settings, channel mapping, and run protocol options.

When the instrument is connected, the Instrument Type and Attune™ Model field are automatically selected based on the instrument and model in use.

Hardware / Virtual Laser Configuration

Instrument Type: Attune NxT Acoustic Focusing Cytometer

Attune Model: 0A28998 Attune NxT Acoustic Focusing Cytometer (Lasers: BVGX)

Autosampler Model: CytKick™ Autosampler

### Instrument type

- **Instrument Type** opens the instrument type dropdown list, which contains the instrument configurations listed in the database.
- Available instrument types are **Attune™ NxT Acoustic Focusing Cytometer** and **Attune™ CytPix™ Flow Cytometer**.

### Attune™ model

- **Attune™ Model** opens the **Instrument configuration dropdown**, which contains the instrument configurations listed in the database.
- The available models are organized by instrument type and the number of lasers in descending order. Within each grouping, models are listed in ascending order.
- A user can import an experiment created on any Attune™ NxT or Attune™ CytPix™ Flow Cytometer model for analysis in the system. However, acquisition is only possible using the model that the system is set to.
- When the system is offline, the **Instrument Configuration dropdown** remains inactive but visible to all user accounts except for the System Administrator and Service accounts, which can modify and set the configuration.
- System Administrator and Service accounts can modify and set the configuration when the Attune™ instrument is offline. This resets the state of the dropdown so that any user can modify the configuration (i.e., if either the System Administrator or Service modifies this, then the dropdown list becomes active for all users).

### Autosampler model

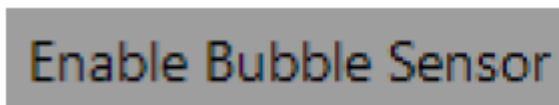
- **Autosampler Model** opens the autosampler model dropdown list, which contains the instrument configurations listed in the database.
- Available autosampler models are **Attune™ Autosampler**, **CytKick™ Autosampler**, and **CytKick™ Max™ Autosampler**.

Autosampler Model:

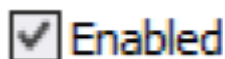
- Attune™ Autosampler
- Attune™ Autosampler
- CytKick™ Autosampler
- CytKick™ Max Autosampler

## Enable bubble sensor

- **Enable Bubble Sensor** option enables the bubble sensor for the system.



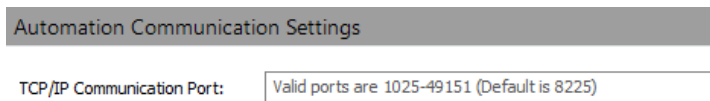
- By default, the bubble sensor option is enabled. This setting is applied to all users.



- Only Administrator, System Administrator, and Service™ accounts can access the Enable Bubble Sensor option. For all other user types, this option is visible but disabled.

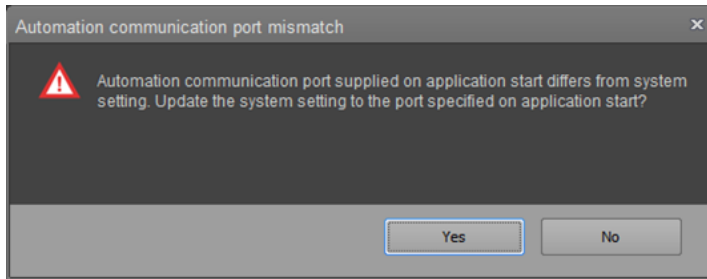
## Automation communication settings

- **TCP/IP Communication Port** allows you to set the port used for communication to and from the automation software (such as the Thermo Scientific™ Momentum™ Software).



- The text box allows values from 1025 to 49151. If the input is outside this range, the text box resets to the last *good* setting, and a warning balloon displays the following text: “Invalid port specified. Valid ports are 1025-49151”.
- To disable the communication port, leave the field blank.
- Click **OK** within the dialog box to set the TCP/IP Communication Port value. Otherwise the changes made to port settings are not preserved.  
Click **Cancel** to revert the port setting to the existing setting. The changes are only applied when the application is restarted.
- Only Administrator, System Administrator, and Service™ accounts can modify the TCP/IP Communication Port settings. For all other user types, this option is visible but disabled.
- When using the Orbitor™ RS3 Microplate Mover with Attune™ NxT and Attune™ CytPix™ instruments, the TCP/IP Communication Port must be set to 8225.
- If the application is started with the port command line argument and the value is different than the application setting, the command line value takes priority.

- When you sign in as an Administrator or System Administrator and there is an Automation communication port mismatch, the *Automation communication port mismatch dialog* allows you to update the application setting.



## Connect device registration

*Connect Device Registration* allows the Administrator or System Administrator to register a device (Attune™ instrument or software) with the Connect cloud-based platform. The Device Name field and the Register Device buttons are only enabled when an Attune™ Administrator or System Administrator is signed into the Attune™ Cytometric Software.

Thermo Fisher Cloud Device Registration

Device Name:

- **Device Name** allows an Administrator or System Administrator to assign a name to the instrument used on the Connect platform when registering the instrument. By default, the device name uses the serial number of the instrument. If this is not available, the machine ID is used as the device name.
- **Register Device** opens the *Link Account* dialog (“Link Account dialog” on page 782). Only Administrators and System Administrators can register a device to the Connect platform. After the device is first registered, different users can log in and log out of their Connect account using the Cloud sign in function (“Sign in to Cloud” on page 53). If the device is already registered, the Register Device button is labeled **Unregister Device**.
- After the instrument has been registered, the Device Name field displays the current device name and the *Update Device Name* button.

Device name was successfully updated.

- To update the device, type in the new name into the Device Name field and click **Update Device Name**. Note that you must have an Administrator or System Administrator account to update the device name.

## Unregister a device from the Connect

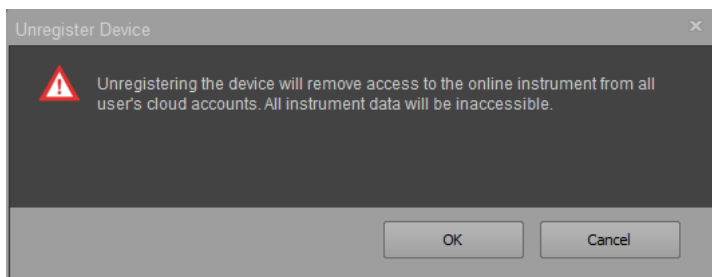
You can completely unregister a device (Attune™ instrument or Attune™ Cytometric software) from the Connect cloud-based platform or dissociate the device from individual Connect accounts.



To completely unregister the device from all Connect accounts, click **Unregister Device**.



The software displays a warning dialog informing that that unregistering the device removes the instrument from the Connect accounts of all users.



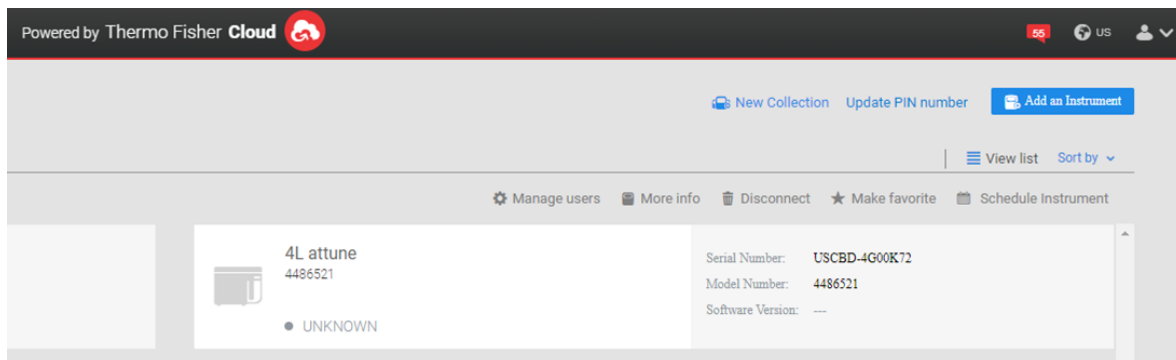
Unregistration completely removes the instrument from the Connect platform as well as from all linked accounts and all data associated with them.

The accounts that were linked to the unregistered Attune™ instrument still exist in the Connect platform and are active for any other devices to which they are connected, but the unregistered Attune™ instrument is no longer listed on the accounts.

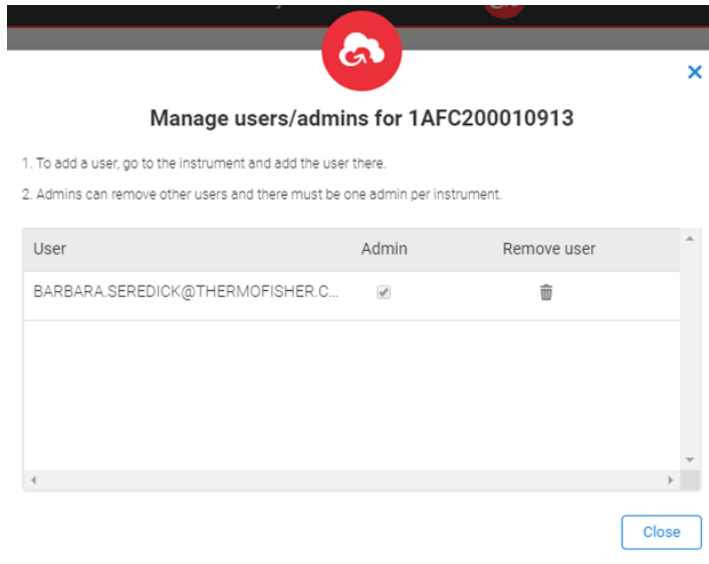
## Remove individual users from a device on the Connect

To remove an individual user from the device, dissociate the device from the Administrator or System Administrator's Connect cloud-based platform account online.

1. Login to the Administrator or System Administrator's Connect account.
2. Within the instrument dashboard, click on the device and select **Manage users**.



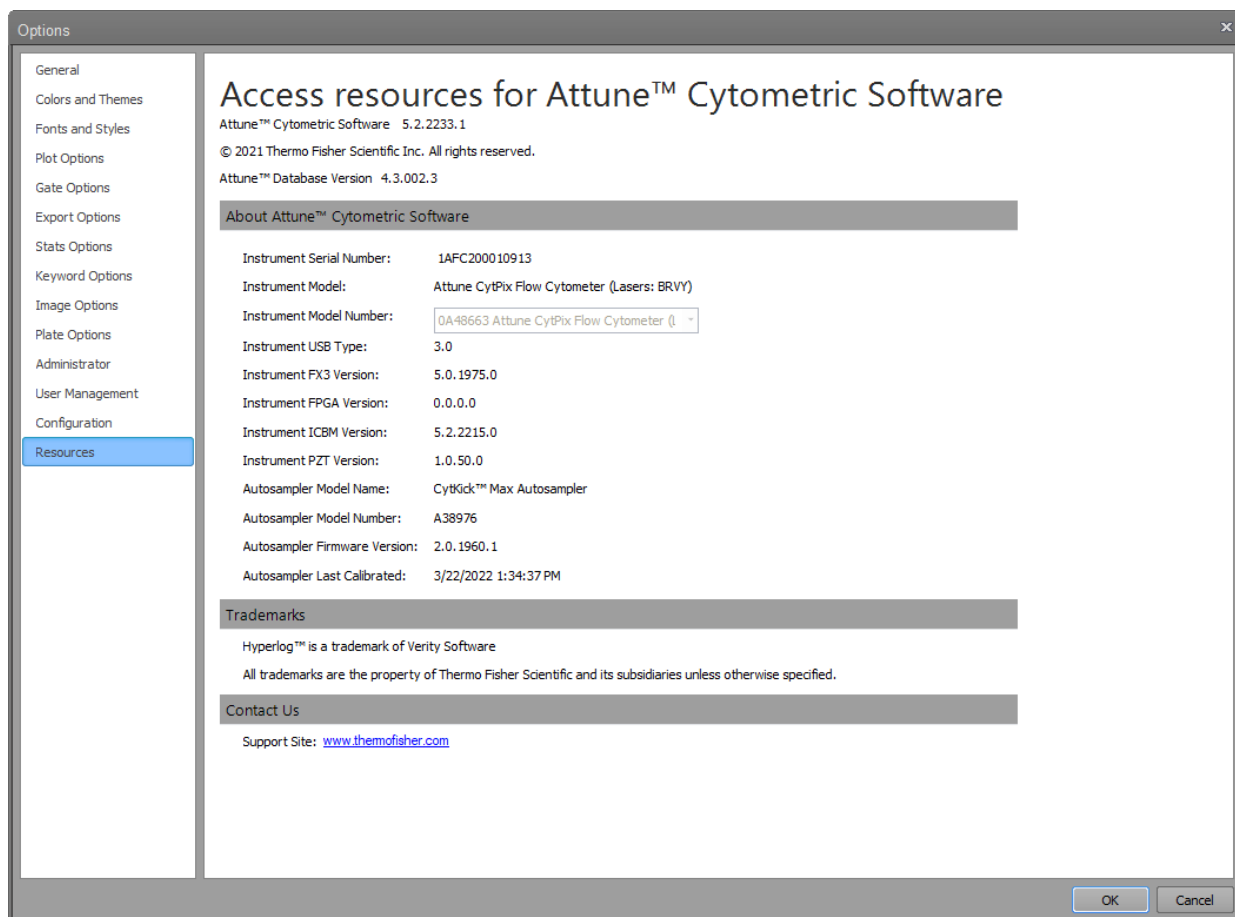
3. In the Manage users/admins dialogue, select the user account to dissociate from the instrument and click on the symbol to remove the user.



# Resources

## Overview

*Resources* tab provides information about the system including the software version, the instrument and auto sampler serial numbers, the instrument and auto sampler firmware versions, the system model, and other licensing and contact information.



## Access resources

This section displays Attune™ Cytometric Software and database version information.

### Attune™ cytometric software

The current Attune™ Cytometric Software version is shown including the software name and software revision number in the format A.B.C.D where A is the major build, B is the minor build, C is the build number and D is the revision.

### Attune™ database version

The current Attune™ database version is read from the database and includes the MySQL version in the format MySQL #.#.#.# | Database #.#.#.#####.

## About Attune™ cytometric software

This section contains the instrument serial number, the instrument model and model number, the instrument firmware, the auto sampler serial number, the auto sampler firmware, and the last calibrated date and time of the auto sampler.

### Instrument serial number

The instrument serial number is automatically retrieved from the instrument.

### Instrument model

- The content for instrument model is automatically populated upon connection to an instrument.
- If no instrument has been connected, the model selected in the *Configuration tab* (“Configuration” on page 695) is shown.
- The instrument model field is updated based on the selection made for instrument model number.

### Instrument model number

- The instrument model number combo box options are automatically populated from the database.
- Upon connection to an instrument, the appropriate model number is automatically selected from the instrument model number combo box options.
- The combo box is disabled by default. It is enabled only for users with Service™ accounts (“Account permissions” on page 44).

### Instrument firmware

- The content for instrument firmware is automatically populated upon connection to an instrument. If no instrument is connected, this field is left blank.
- The instrument firmware field displays the major, minor, and build revision number in the format `##.##.###.#`.

### Auto sampler serial number

- The auto sampler serial number field is blank by default. A user with a Service™ account can enter an auto sampler serial number.
- The maximum length for a serial number is 50 characters. Any ASCII characters can be entered except `/ \ ? % * : | ” < >`.  
Spaces are not permitted and cannot be entered.
- If a user attempts to enter an illegal character, a warning dialog indicates the error condition, and the invalid characters are not entered.

### Auto sampler last calibrated

- The section is automatically updated to show the date and time of the last auto sampler calibration. This value is read from the instrument.
- The data and time are displayed in the local date and time format.

## Trademarks

This section contains information on any trademarks used within the software.

## Contact us

This section contains hyperlinks to Thermo Fisher Scientific websites.

**User Management** tab of the **Options dialog** enables authorized users to perform the following user management functions:

- Add or edit a local Attune™ account (“Add or edit an account” on page 705)
- Reset a password for a local Attune™ account (“Reset a password” on page 711)
- Change a password for a local Attune™ account (“Change a password” on page 712)

---

**Note:** When the **SAE mode** is enabled (see “Enable the SAE mode” on page 682), user management is controlled by the SAE Administrator Console (see Appendix D, “SAE Administrator Console”). For SAE specific permissions, see “SAE account permissions” on page 887.

---

## Access the user management tab

1. In the **Quick Access** toolbar, click **Options** to open the **Options dialog**.
2. In the left pane, select the **User Management** tab.

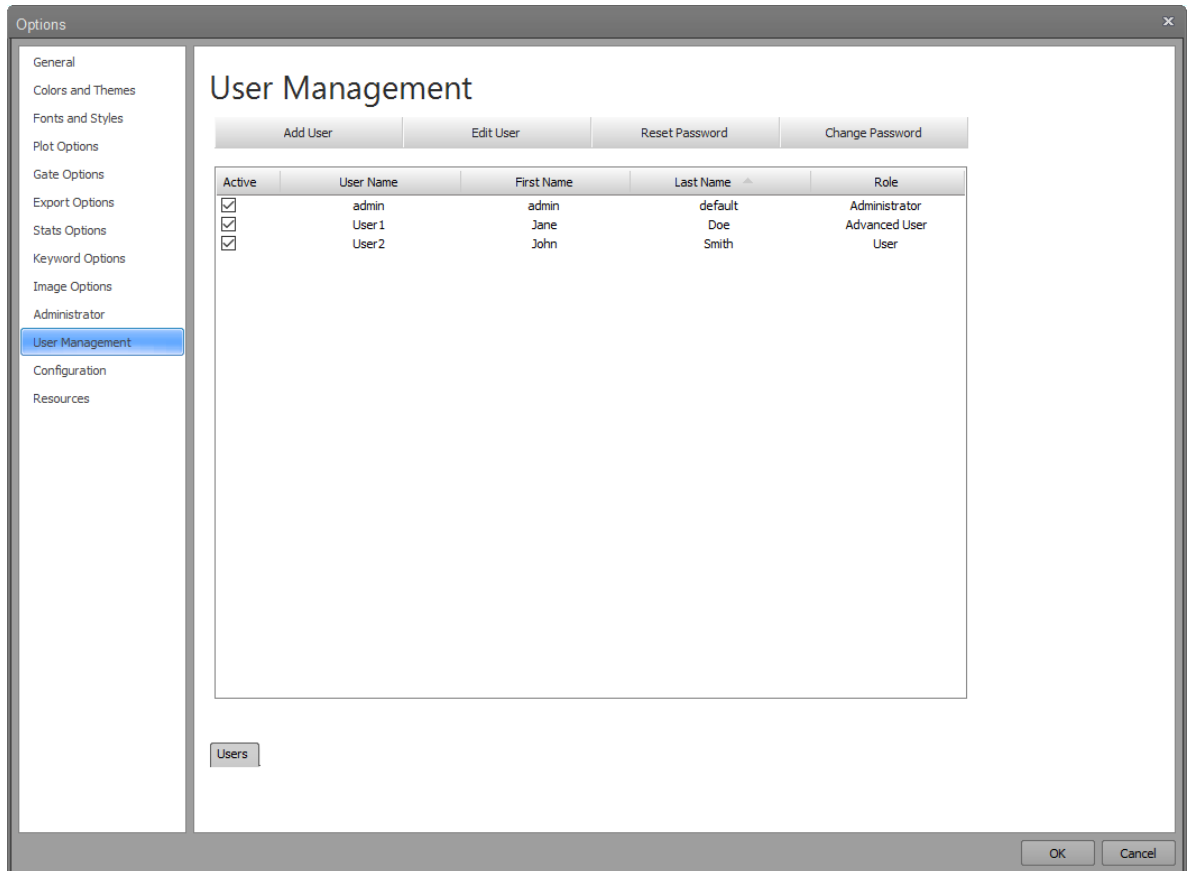


Figure 137 User Management tab – Administrator view

## Add or edit an account

### Add an account

Only **System Administrator** and **Administrator** accounts (“Default local Attune™ accounts” on page 42) can add new local Attune™ accounts to the software. To create a user profile to add a new account, see “Create a user profile” on page 706.

### Edit an account

**System Administrator** and **Administrator** accounts can edit all local Attune™ accounts (except the **Service** account, which cannot be edited). **Users** can only edit their own accounts. To edit an existing user account, see “Edit a user profile” on page 708.

## Create a user profile

Only **System Administrator** and **Administrator** accounts can create a local Attune™ user profile.

1. In the **User Management** tab of the **Options dialog**, click **Add User** to open the **Add User dialog**.

**Add User** [Close]

**Profile**

User Name\*:

First Name\*:

Last Name\*:

Phone Number:

Email Address:

Principal Investigator:

Select Roles:

Permissions Bases on Selected Roles

- Delete Report Column Access
- Adjust Laser Delays
- Database Utility Full Access
- View Shared Folders
- Run Performance Test
- Advanced Instrument Settings
- Service User Tools
- Run System Decontamination
- System Instrument Settings
- Manage User Accounts
- Modify Instrument Configuration
- Run System Tests

OK Cancel Help



2. Enter the requested profile information:

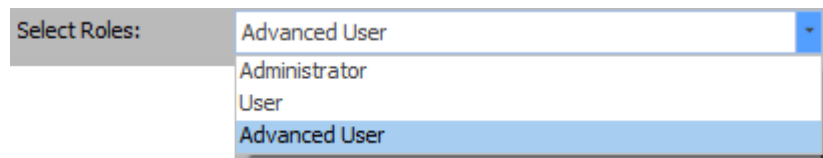
- **Username\***: Required field. Enter a unique username between 3 and 50 characters. Duplicate usernames are not permitted. Spaces and the following characters are not permitted: \ / : \* ? " < > |
- **First Name\***: Required field. Enter up to 50 characters; the following characters are not permitted: \ / : \* ? " < > |
- **Last Name\***: Required field. Enter up to 50 characters; the following characters are not permitted: \ / : \* ? " < > | %
- **Phone Number**: Optional field. Enter up to 50 characters. The following characters are permitted: +, and the numerals 0 through 9  
The + character can only be in the first position; for example: **+0123456789**
- **Email Address**: Optional field. Enter up to 100 characters; the following characters are not permitted: \ / : \* ? " < > | %  
The email address must contain the @ character and a domain name; for example, **email@address.com**.
- **Principal Investigator**: Optional field. Enter up to 100 characters; the following characters are not permitted: \ / : \* ? " < > | %

---

**Note:** For all fields, the software automatically removes trailing, leading, or consecutive spaces.

---

3. Select an account type from the **Select Roles** dropdown menu: **User**, **Advanced User**, **Administrator**.



- You cannot add a new **Service** account to the system.
  - Permissions assigned to the selected account are displayed in black text; permissions not assigned to the account are grey. None of the user account permissions can be modified.
  - For a list of account permissions, see “Default local Attune™ accounts” on page 42.
4. Click **OK** to create the user account and close the dialog.  
To close the dialog without creating an account, click **Cancel** or **X**.

---

**Note:** For the permissions assigned to each local Attune™ account type, see “Account permissions” on page 44.

For SAE specific permissions, see “SAE account permissions” on page 887.

---

## Edit a user profile

**System Administrator** or **Administrator** accounts can edit all local Attune™ accounts (except the **Service** account, which cannot be edited). **Users** can only edit their own local Attune™ accounts. **SAE users** cannot edit their user profiles using the Attune™ Cytometric Software and must use the SAE Administrator Console (Appendix D, “SAE Administrator Console”).

**Note:** For the permissions assigned to each local Attune™ account type, see “Account permissions” on page 44.

For SAE specific permissions, see “SAE account permissions” on page 887.

1. In the **User Management** tab of the **Options dialog** (“Users tab” on page 688), click **Edit User** to open the **Edit User dialog**.

**Edit User**

**Profile**

User Name\*: User 1

First Name\*: Jane

Last Name\*: Doe

Phone Number:

Email Address:

Principal Investigator:

Select Roles: Advanced User

Permissions Bases on Selected Roles

Delete Report Column Access	Service User Tools
Adjust Laser Delays	Run System Decontamination
Database Utility Full Access	System Instrument Settings
View Shared Folders	Manage User Accounts
Run Performance Test	Modify Instrument Configuration
Advanced Instrument Settings	Run System Tests

OK Cancel Help

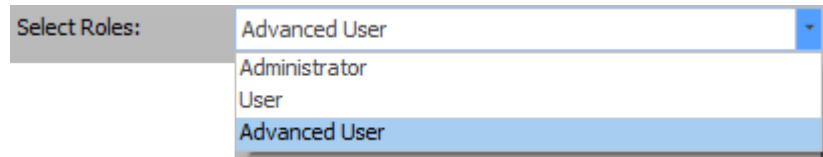
2. In the **Profile** tab, edit the desired profile information fields as described in “Add or edit an account” on page 705.

---

**Note:** For all fields, the software automatically removes trailing, leading, or consecutive spaces.

---

3. Select an account type from the **Select Roles** dropdown menu: **User**, **Advanced User**, or **Administrator**.

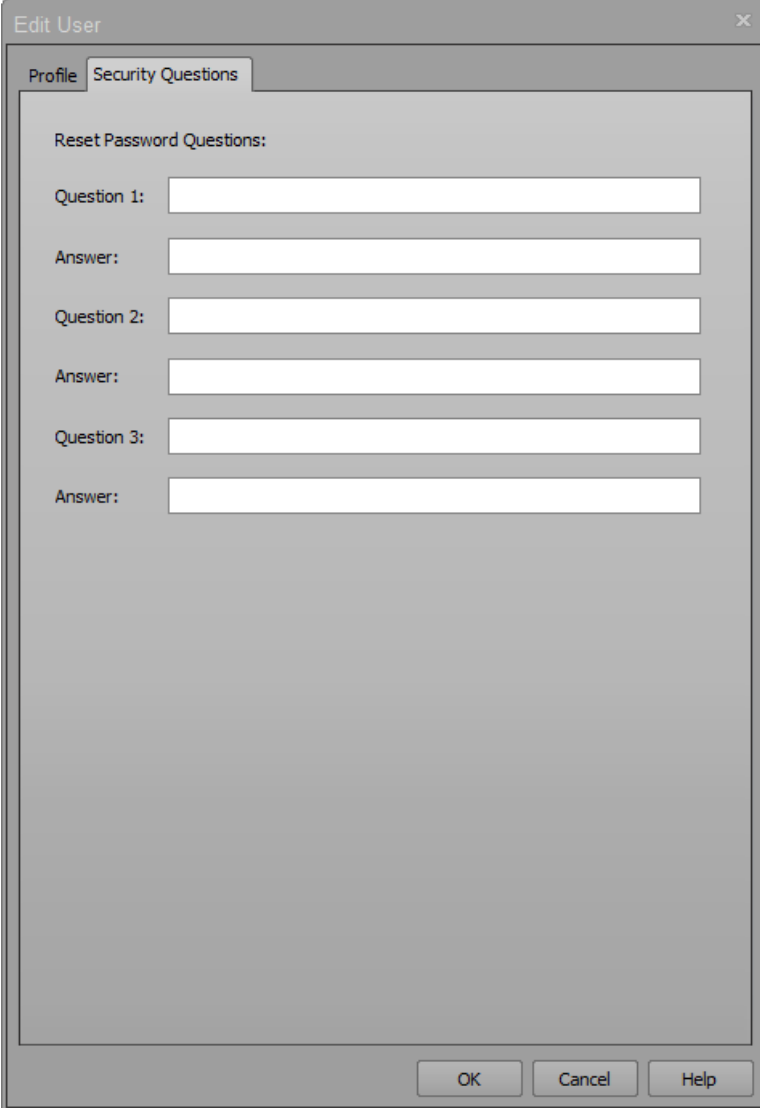


- Only **System Administrator** or **Administrator** accounts can edit the account type. The **Select Roles** dropdown is disabled for all other users.
  - Permissions assigned to the selected account are displayed in black text; permissions not assigned to the account are grey. None of the user account permissions can be modified.
  - For a list of user account permissions, see “Default local Attune™ accounts” on page 42.
4. Click **OK** to save the changes and close the dialog.  
To close the dialog without saving the changes, click **Cancel** or **X**.

## Set password reminders

You can set password reminders only for your own local Attune™ account.

1. In the **User Management** tab, click **Edit User** to open the **Edit User dialog**.
2. In the **Edit User dialog**, click the **Security Questions** tab.



The screenshot shows the 'Edit User' dialog box with the 'Security Questions' tab selected. The dialog has a title bar with 'Edit User' and a close button. Below the title bar are two tabs: 'Profile' and 'Security Questions'. The 'Security Questions' tab is active and contains the following fields:

Reset Password Questions:

Question 1:

Answer:

Question 2:

Answer:

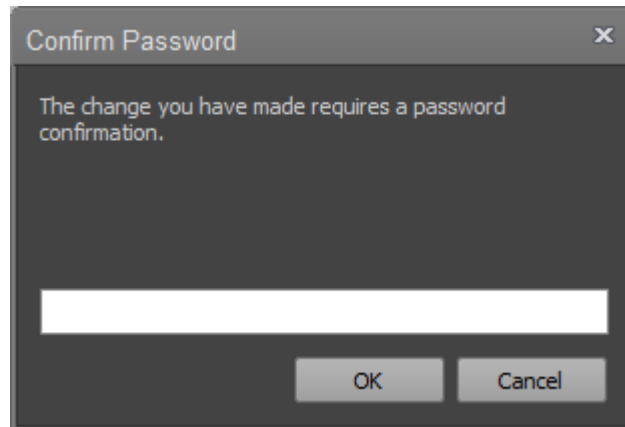
Question 3:

Answer:

At the bottom of the dialog are three buttons: 'OK', 'Cancel', and 'Help'.

3. Enter up to three password reminder questions and answers, then click **OK**.
  - For each question and answer, you can enter up to 100 characters; any character is permitted.
  - To prevent unauthorized access to a user account, the answers are displayed as bullet points.
  - You can reset a forgotten password by correctly answering the password reminder questions.

- At the prompt, enter your password to confirm the password reminder changes, then click **OK**.

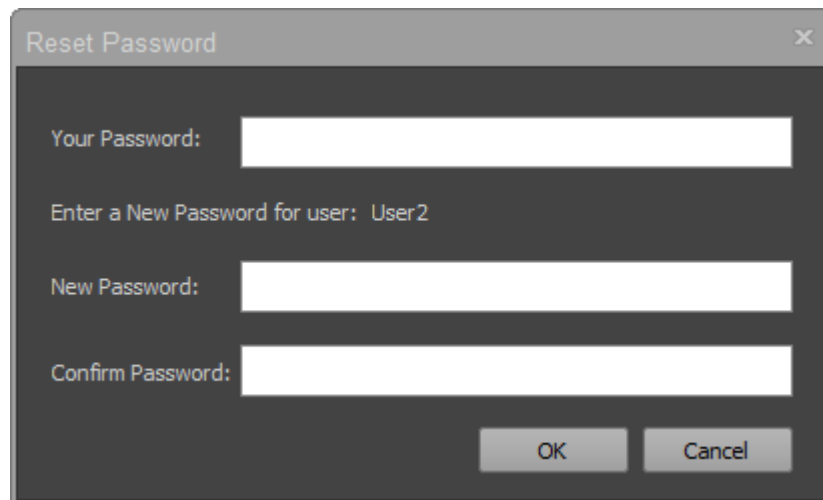


- In the **Edit User dialog**, click **OK** to save the changes and close the dialog. To close the dialog without saving the changes, click **Cancel** or **X**.

## Reset a password

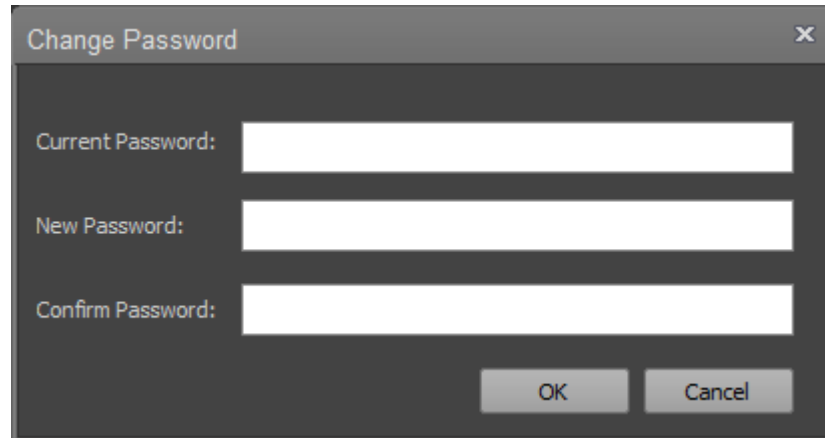
**Reset Password** provides **System Administrators** and **Administrators** the ability to reset the password for local **User** and **Advanced User** accounts. Only **System Administrator** or **Administrator** accounts can reset a password. **User** and **Advanced User** accounts cannot reset the passwords of other accounts.

- In the **User Management** tab, click **Reset Password** to open the **Reset Password dialog**.



- For **Your Password**, enter the System Administrator or Administrator password.
- For **New Password**, enter the new password for the account whose password is to be reset.
- For **Confirm Password**, re-enter the new password.

5. Click **OK** to save the changes and close the dialog.  
To close the dialog without saving the changes, click **Cancel** or **X**.
6. Provide the user with their new password. At next login, the user will be prompted to change the password.



---

**Note:** For local Attune™ users, the password policy applied to reset passwords is based on the system security policy for user passwords set in the **User Management ▶ System Security** tab of the **Options dialog** (see “System Security tab” on page 692).

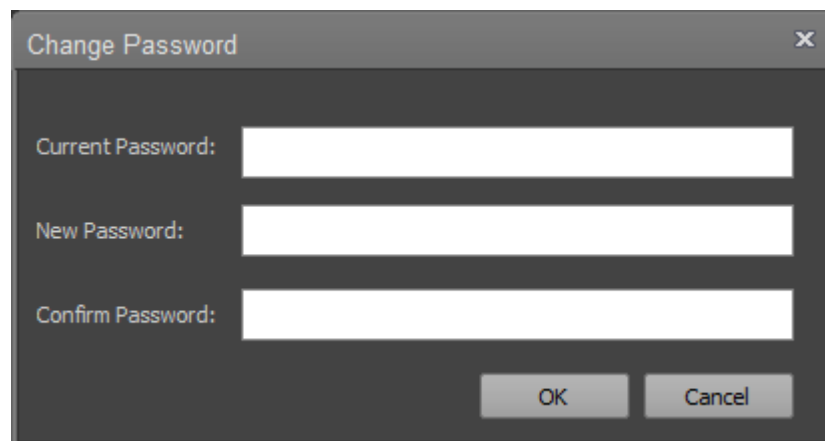
---

For SAE users, the password policy applied to reset passwords is set based on the SAE security policy (see “SAE account permissions” on page 887).

## Change a password

**Change Password** provides any user the option to change their own local Attune™ password.

1. In the **User Management** tab, click **Change Password** to open the **Change Password dialog**.



2. For **Current Password**, enter your existing password.

3. For **New Password**, enter a new password, then confirm the new password in **Confirm Password** field.
4. Click **OK** to save the changes and close the dialog.  
To close the dialog without saving the changes, click **Cancel** or **X**.

---

**Note:** For local Attune™ users, the password policy applied to changing passwords is based on the system security policy for user passwords set in the **User Management ▶ System Security** tab of the **Options dialog** (see “System Security tab” on page 692).

For SAE users, the password policy applied to changing passwords is set based on the SAE security policy (see “SAE account permissions” on page 887).

---

## Overview

This chapter describes the dialogs used for opening and saving files and folders, loading configuration files (Workspace, Instrument Settings, Heat Map, Run Protocols, Compensation Settings) stored in the database, and dialogs that are displayed when importing or exporting FCS files and images.

All dialogs are modal, unless otherwise specified.

---

**Note:** Modal dialogs appear when you are performing tasks in the main software windows. You must interact with the modal dialog before you can return to the main window.

---

## File dialogs

- **File Save (Export)** dialog (“File Save (Export) dialog” on page 715)
- **File Open (Import)** dialog (“File Open (Import) dialog” on page 721)
- **Folder Browser** dialog (“Folder Browser dialog” on page 723)



# File Save (Export) dialog

## Overview

**File Save (Export)** dialog uses the standard Microsoft™ Windows™ browser, which enables you to save the selected files in the desired location. Depending on the file that is being saved (for example Experiment, FCS file, image file, sample list, etc.), the dialog provides different save options (for example FCS file version, image format, etc.).

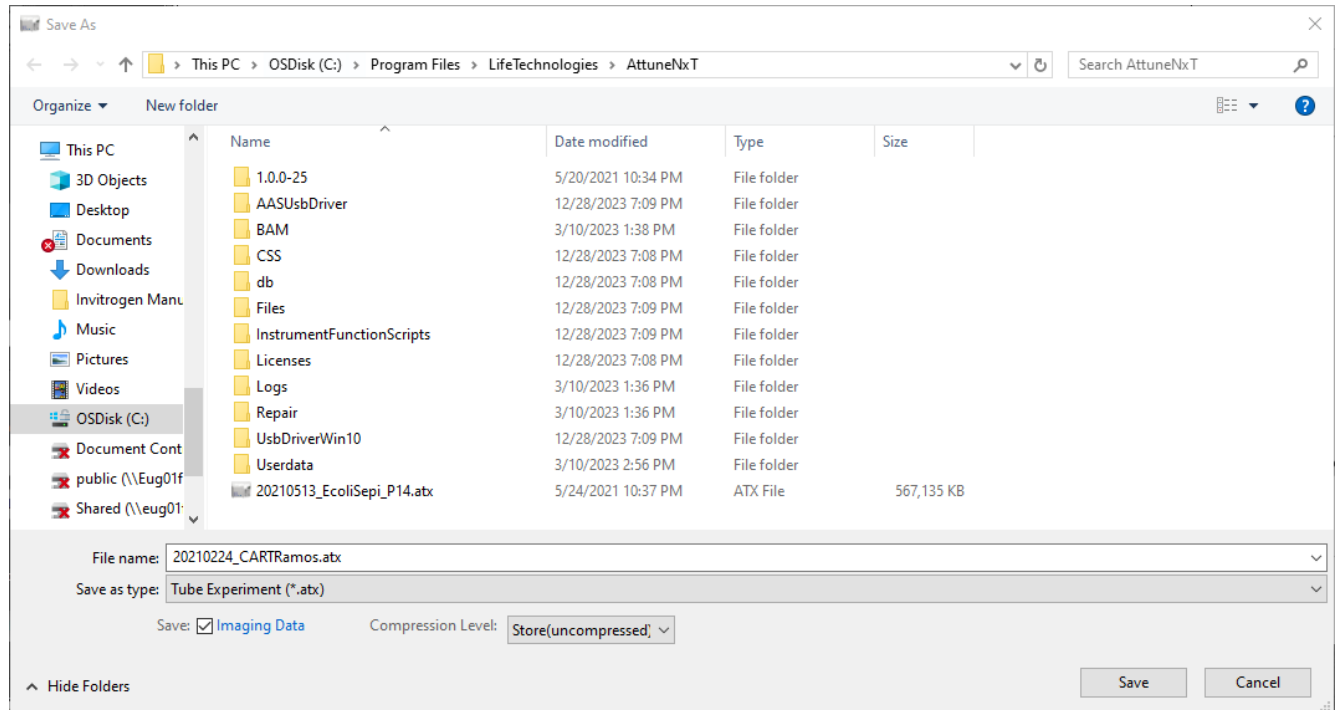
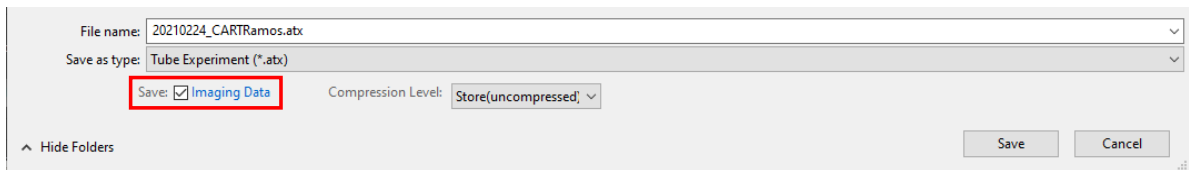


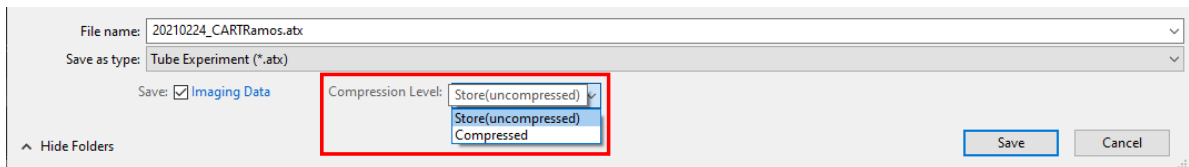
Figure 138 File Save (Export) dialog for saving a Tube Experiment with the option to save imaging data.

## Dialog behavior

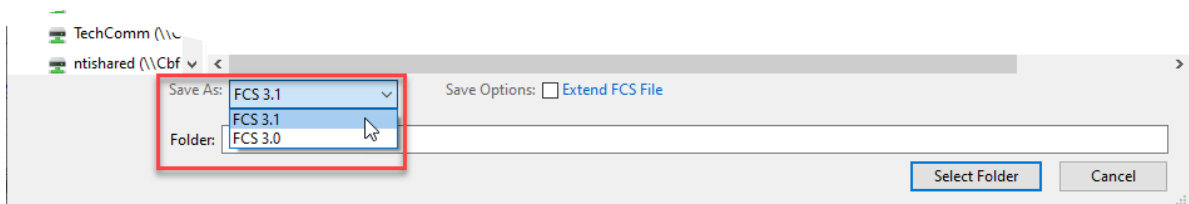
- The **File Save (Export)** dialog opens to the last visited folder where files were saved.
- Depending on the file that is being saved (for example Experiment, FCS file, image file, sample list etc.), the dialog provides different save options (for example FCS file version, image format, compression level etc.).
- When saving, the file extension is automatically added to each file type and the **Confirm File Save** dialog (“Confirm file save dialog” on page 746) is displayed, if needed.
- If a name with invalid characters is entered or a name that results in a path length over 260 characters, a warning dialog indicating the error condition is displayed.
- Clicking **OK** on the warning dialog returns to the **File Save (Export)** dialog and the file is not saved.
- When exporting an Experiment that has captured images, you can save the imaging data together with the Experiment metadata and sample data by selecting the **Save Imaging Data** option.



- When exporting an Experiment or images, you can save the data files uncompressed by selecting **Store (uncompressed)** from the **Compression Level** dropdown. Alternatively, you can select **Compressed** to save storage space when exporting the Experiment or the images.

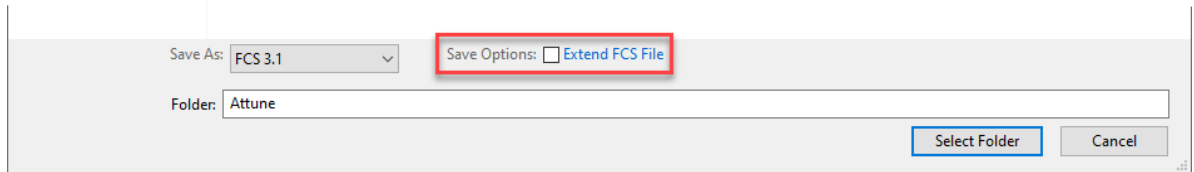


- When saving FCS files, the default filter is set to the **FCS 3.1** version. You can save files as FCS 3.0 by selecting **FCS 3.0** from the **Save As** dropdown (when exporting from an Experiment or Group) or from the **Save as type** dropdown (when exporting from a Sample).



- When exporting sample data that has image processing data, the image processing results can be appended to the original FCS file, which creates a single FCS file that includes the image processing parameters.

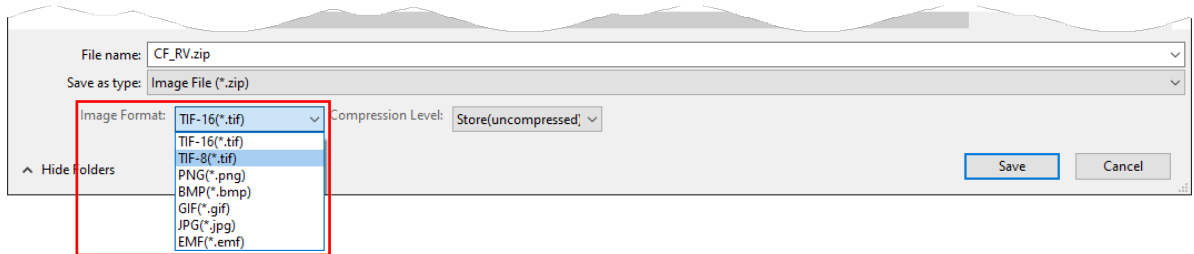
To export FCS files and imaging data together, select **Extend FCS File** for **Save Options** to add the new data to the original FCS file. This option can be used for whole datasets or specifically for exporting from a gate.



When **Extend FCS File** option is selected, the original FCS file is appended with the image processing parameter data, and the \$ORIGINALITY, \$LAST\_MODIFIER, and \$LAST\_MODIFIED keywords in the FCS file are updated to indicate that the data have been appended and modified.

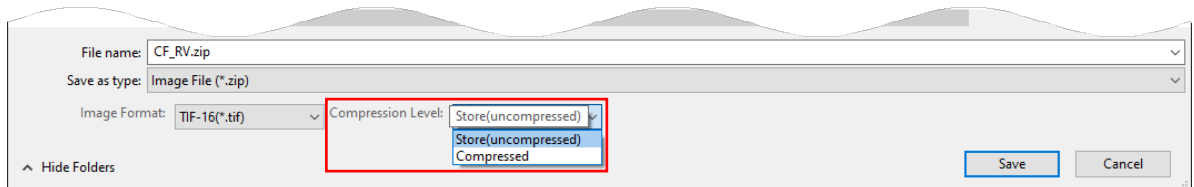
- When **Export Images** is selected to archive captured images from an Experiment, Group, or Sample, all images are exported into a single zip file.

By default, images are exported as **TIF-16** files, which is the native file format collected from the Attune™ CytPix™ Flow Cytometer. You can select to save the images as **TIF-16**, **TIF-8**, **PNG**, **BMP**, **GIF**, **EMF**, or **JPG** files using the **Image Format** dropdown.



**IMPORTANT!** When saving in any format other than TIF-16, images are converted to 8-bit images, which results in loss of image pixel data.

The **Compression Level** dropdown enables you to export the image files as **Store (uncompressed)** or as **Compressed**.



## File filters and default names

The file filters and default names depend on the source of the File Save (Export) dialog and are described in the following table.

Action/Source	Default Extension	Default File Name	Filter
Export FCS File	FCS	Name of File as saved by user	FCS 3.0 (*.fcs) FCS 3.1 (*.fcs)
Export FCS and Image Data	ACS	Name of File as saved by user	Archival Cytometry Standard (*.acs)
Export Heat Map	AHM	Plate_Experiment Experiment	HeatMap File (*.ahm)
Export Run Protocol	ARP	Plate_Experiment_Group_Sample Experiment_Group_Sample	Run Protocol File (*.arp)
Export Workspace	AWS	Experiment-level: Plate_Experiment Group-level: Plate_Experiment_Group Sample-level: Plate_Experiment_Group_Sample	Workspace File (*.aws)
Export Instrument Settings	AIS	Experiment-level: Plate_Experiment Sample-level: Plate_Experiment_Group_Sample	Instrument Settings File (*.ais)
Export Template	AET	Tube or Plate	Template (*.aet)
Export Plate Experiment	APX	Plate	Plate Experiment (*.apx)
Export Tube Experiment	ATX	Experiment	Tube Experiment (*.atx)
Export Statistics (Single File Output)	CSV	Sample-level: Plate_Experiment_Group_Sample Results tab: Plate_Experiment Multiple files: "Blank"	CSV (*.csv)
Export Compensation Settings	ACS	Experiment-level: Experiment Sample-level: Plate_Experiment_Group_Sample	Compensation Settings (*.acs)
Send Table to XML	XML	Name of file as saved by user	XML (*.xml)

(continued)

Action/Source	Default Extension	Default File Name	Filter
Print to PDF	PDF	Sample-level: Plate_Experiment_Group_Sample Results tab: Plate_Experiment Multiple files: "Blank"	PDF (*.pdf)
Save as Powerpoint	PPTX	Report: Plate_Experiment_Group_Sample	Powerpoint (*.pptx)
Save System Log	ALF	Long Date in system locale format	Log File (*.af)
Export Gate to FCS	FCS	Sample_GateName	FCS 3.0 (*.fcs) FCS 3.1 (*.fcs)
Export Images	ZIP	Name of File as saved by user	TIF-16 (*.tif) TIF-8 (*.tif) PNG (*.png) BMP (*.bmp) GIF (*.gif) JPG (*.jpg) EMF (*.emf)
Save as (to image)	PNG	Sample	PNG (*.png) GIF (*.gif) JPG (*.jpg) TIF (*.tif) BMP (*.bmp) PDF (*.pdf) Windows™ Metafile
Save as (cell image)	TIF-8	"Blank"	TIF-8 (*.tif) TIF-16 (*.tif) PNG (*.png) BMP (*.bmp) GIF (*.gif) JPG (*.jpg) EMF (*.emf)
Export Instrument Configuration	AIC	Instrument Configuration Name	AIC (*.aic)

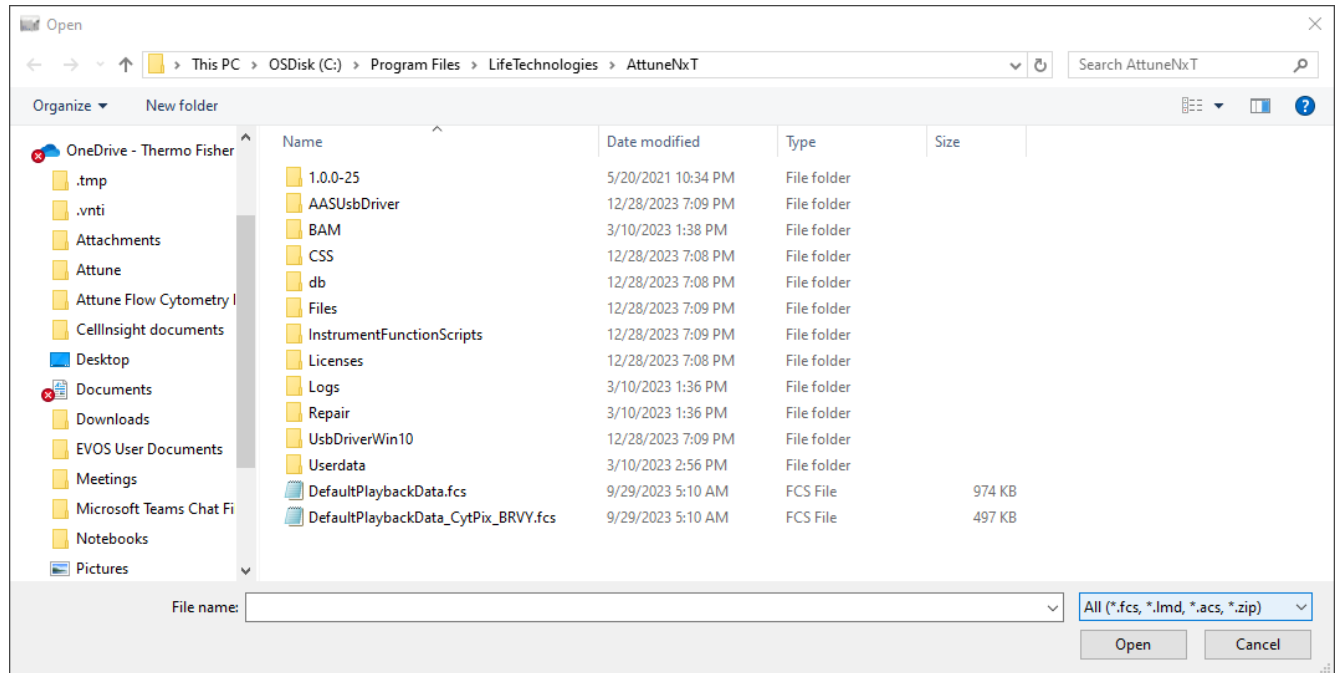
(continued)

Action/Source	Default Extension	Default File Name	Filter
Export Plates	APL	Multi-select: "Attune™ Plate Definitions" Single-select: <plate name>	APL (*.apl)
Sample List	CSV	Plate or Experiment name	CSV (*.csv)
Export Logs for Service	ZIP	AttuneServiceLogs_YYYY-MM-DD_HH-MM-SS (where YYYYMMDD_HHMMSS is the date and time)	ZIP (*.zip)

# File Open (Import) dialog

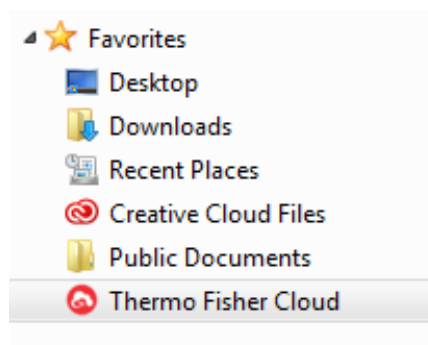
## Overview

**File Open (Import)** dialog uses the standard Microsoft™ Windows™ File Open browser. It is displayed when **Import** is selected in the software.



## Dialog behavior

- The **File Open (Import)** dialog opens to the last visited folder with open files.
- FCS files, Plate and Tube Experiments can be imported.
- If the device (instrument or software) has been registered to a Connect account, the **File Open (Import)** dialogue includes an icon for the Connect listed under **Favorites**. Clicking on this icon opens a virtual folder listing all files and folders included in the Connect account.



- Multiselect is enabled for FCS files when **Import FCS Files** is selected from the **Experiment context menu** on Experiment Explorer (“Rename Experiment” on page 312), using the New Experiment dialog’s Analysis type experiment or Compensation Setup Dialog’s “Files” source option. Multiselect is also available for Plate and Tube experiments. Multiselect is not enabled for the other file types.
- When FCS files are imported, more options to update the experiment voltages and experiment compensation are provided.
- If the Instrument settings of the imported FCS file do not match the current Instrument configuration, no option to update the Experiment voltages or experiment compensation is provided.
- When multiple FCS files are selected, no option to update the experiment voltages or experiment compensation is provided

## File filters

The file filters depend on the source of the **File Open (Import)** dialog command and are described in the following table.

Action	Filter
Import FCS File	All (*.fcs,*.lmd)
Import Heat Map	Heat Map File (*.ahm)
Import Run Protocol	Run Protocol File (*.arp)
Import Workspace	Workspace File (*.aws)
Import Instrument Settings	Instrument Settings File (*.ais)
Import Template	Plate Template (*.aet)
Import Experiment files	All (*.apx, *.atx) <b>Note:</b> Plate Experiment (*.apx), Tube Experiment (*.atx)
Import Compensation Settings	Compensation Settings File (*.acs)
Import Bead Lot File	Bead Lot File (*.abl)
Import Instrument Configuration	Instrument Configuration (*.aic)

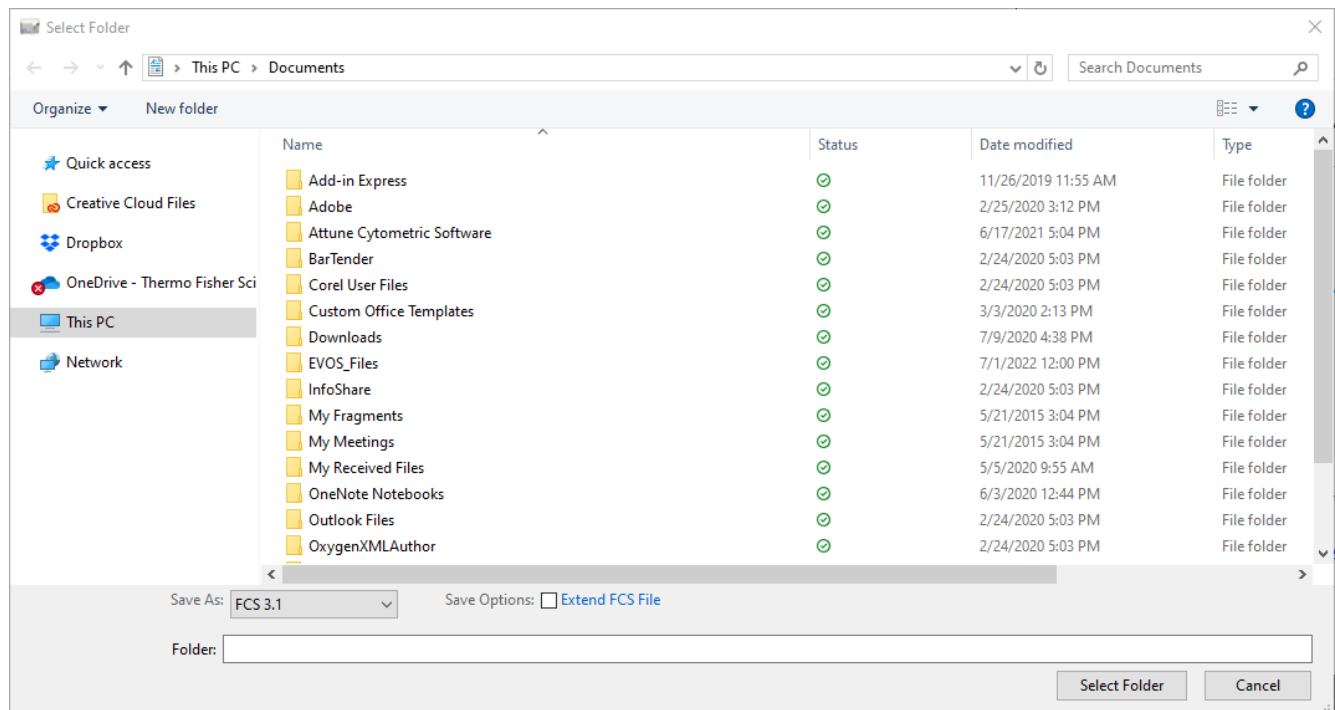


# Folder Browser dialog

## Overview

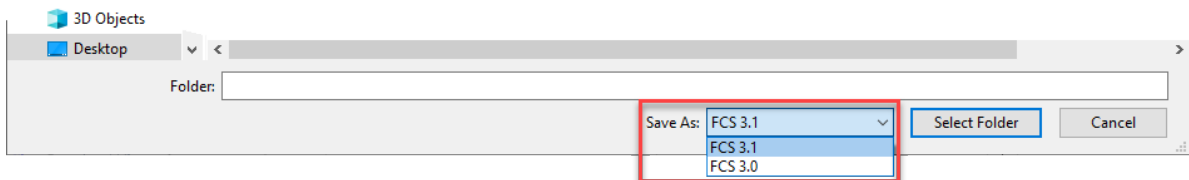
**Folder Browser** dialog is displayed when multiple files can be opened by selecting a folder or multiselecting several files in a folder. Additionally, this dialog is used when a **Save** command requires the selection of a folder.

The style of the **Folder Browser** is the same as those used for the **File Save (Export)** and **File Open (Import)** browsers.

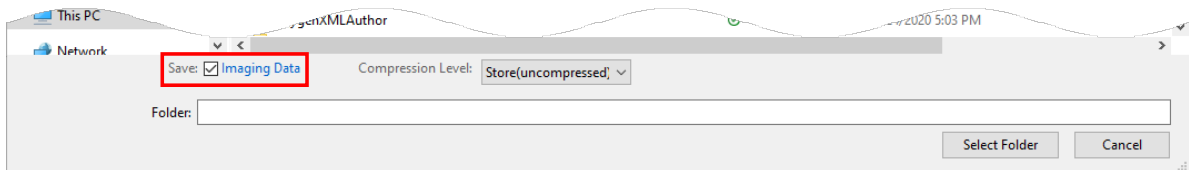


## Dialog behavior

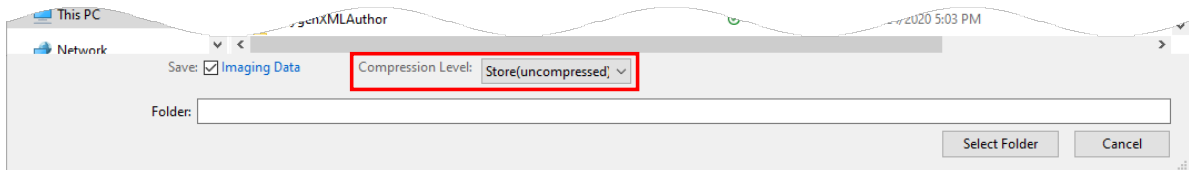
- The **Folder Browser** dialog opens to the last visited folder where files were open.
- Depending on the file that is being saved (for example Experiment, FCS file, image file, sample list etc.), the dialog provides different save options (for example FCS file version, image format, compression level etc.).
- When exporting, the file extension is automatically added to each file type and the **Confirm File Save** dialog (“Confirm file save dialog” on page 746) is displayed for each save action.
- When exporting FCS files, the folder dialog includes the **Save As** option, which lets you export FCS files as **FCS 3.1** or **FCS 3.0** files. The default choice is set to the **FCS 3.1**.



- When exporting multiple Experiments that have captured images, you can save the imaging data together with the Experiment metadata and sample data by selecting the **Save Imaging Data** option.

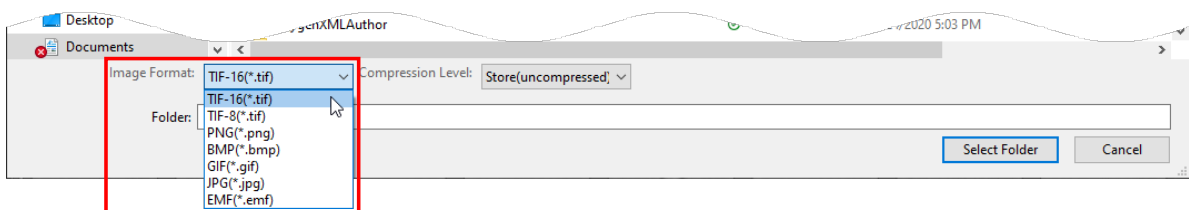


- When exporting an Experiment or images, you can save the data files uncompressed by selecting **Store (uncompressed)** from the **Compression Level** dropdown. Alternatively, you can select **Compressed** to save disk storage space when exporting the Experiment or the images.



- When **Export Images** is selected to archive captured images from multiple Experiments, Groups, or Samples, the images from each Experiment, Group, or Sample are exported into separate zip files that include all images from that Experiment, Group, or Sample.

By default, images are exported as **TIF-16** files, which is the native file format collected from the Attune™ CytPix™ Flow Cytometer. You can select to save the images as **TIF-16**, **TIF-8**, **PNG**, **BMP**, **GIF**, **EMF**, or **JPG** files using the **Image Format** dropdown.

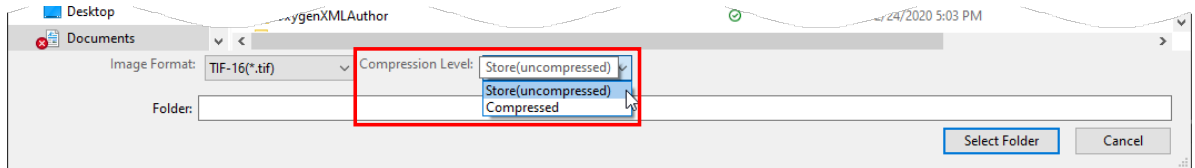


---

**IMPORTANT!** When saving in any format other than **TIF-16**, images are converted to 8-bit images, which results in loss of image pixel data.

---

By default, **Store (uncompressed)** is selected for **Compression Level** when saving images from multiple Experiments, Groups, or Samples. To save disk storage space, you can select **Compressed** from the **Compression Level** dropdown list.



## File filters and default names

When exporting, file filters and default names are as described in the following table.

Action/Source	Default Extension	Default File Name	Filter
Import FCS Files	FCS	Name of File as saved by user	All (*.fcs,*.lmd)
Export FCS Files	FCS	Plate_Experiment_Group_Sample	FCS 3.1 (*.fcs)
Export FCS and Image Data	ACS	Name of File as saved by user	Archival Cytometry Standard (*.acs)
Export Statistics (Multiple File Output)	CSV	Sample-level: Plate_Experiment_Group_Sample	CSV (*.csv)
Export Images	ZIP	Name of File as saved by user	TIF-16 (*.tif) TIF-8 (*.tif) PNG (*.png) BMP (*.bmp) GIF (*.gif) JPG (*.jpg) EMF (*.emf)

## Overview

This chapter describes the non-instrument dialogs that are displayed throughout the Attune™ Cytometric Software.

All dialogs are *modal*, unless otherwise specified.

---

**Note:** Modal dialogs appear when you are performing tasks in the main software windows. You must interact with the modal dialog before you can return to the main window.

---

- Page Setup dialog (“Page Setup dialog” on page 728)
- Print dialog (“Print dialog” on page 730)
- Print Preview dialog (“Print Preview dialog” on page 732)
- Batch Print dialog – Experiment Explorer context menu (“Batch Print dialog” on page 734)
- Low Disk Space dialog (“Low Disk Space dialog” on page 736)
- Export Statistics dialog – Experiment Explorer context menu (“Export Statistics dialog” on page 737)
- Derived Gate dialog (“Derived gate dialog” on page 739)
- Edit Gates dialog (“Edit gates dialog” on page 742)
- Confirm File Save dialog (“Confirm file save dialog” on page 746)
- Save as Template dialog (“Save As Template dialog” on page 747)
- Manage Templates dialog (“Manage templates dialog” on page 749)
- Plate Mapping dialog (“Plate Mapping dialog” on page 756)
- Select Template dialog (“Select template dialog” on page 760)
- Map Sample List Data dialog (“Map Sample List Data dialog” on page 764)
- Update Sample Information dialog (“Update sample information dialog” on page 766)
- Create/Edit Keyword dialog (“Create/Edit keyword dialog” on page 767)
- Experiment Keywords dialog (“Experiment Keywords dialog” on page 770)
- Export Logs for Service dialog (“Export Logs for Service dialog” on page 772)
- Select Plates dialog (“Select Plates dialog” on page 774)
- Create/Edit Plate dialog (“Create/Edit plate” on page 776)
- Test Plate dialog (“Test Plate dialog” on page 778)
- Foil cover present dialog (“Foil cover present dialog” on page 779)
- Process (Reprocess) Images dialog (“Process (Reprocess) Images dialog” on page 780)
- Link Account dialog (“Link Account dialog” on page 782)
- Cloud Sign in dialog (“Cloud Sign In dialog” on page 787)

- Switch Account dialog (“Switch Account dialog” on page 790)
- Connect Storage dialog (“Connect storage dialog” on page 791)
- Maintenance Required dialog (“Maintenance Required dialog” on page 792)
- System Log dialog (“System Log dialog” on page 793)
- SAE signing dialogs (“SAE Signing dialogs” on page 804)

## Page Setup dialog

The **Page Setup dialog** enables you to customize the paper size and orientation.

### Open the dialog

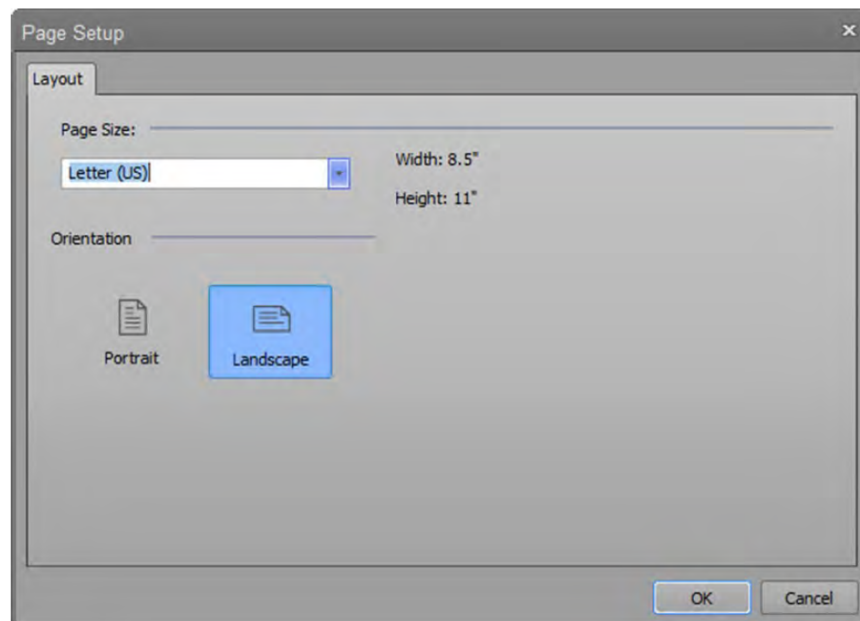
On the **Home ribbon tab** (“Home tab” on page 74), click **Page Setup**.

The **Page Setup** button is only available when the Workspace, Results, Heat Map, or Overlay views are active.

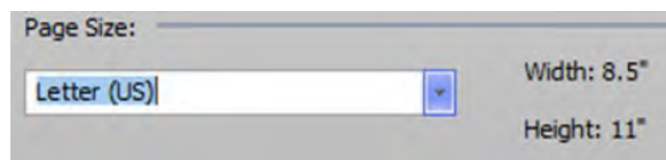


### Layout tab

The **Layout tab** of the **Page Setup dialog** enables you to customize the page size and orientation.

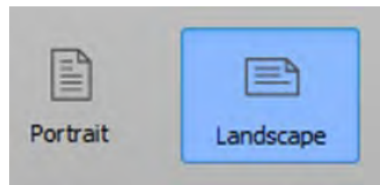


1. To select a page size for the printed document, select a size from the **Page Size** dropdown menu.



The software displays the width and height dimensions for the selected size.

2. To select the orientation of the printed document, click the **Portrait** or **Landscape** icon.



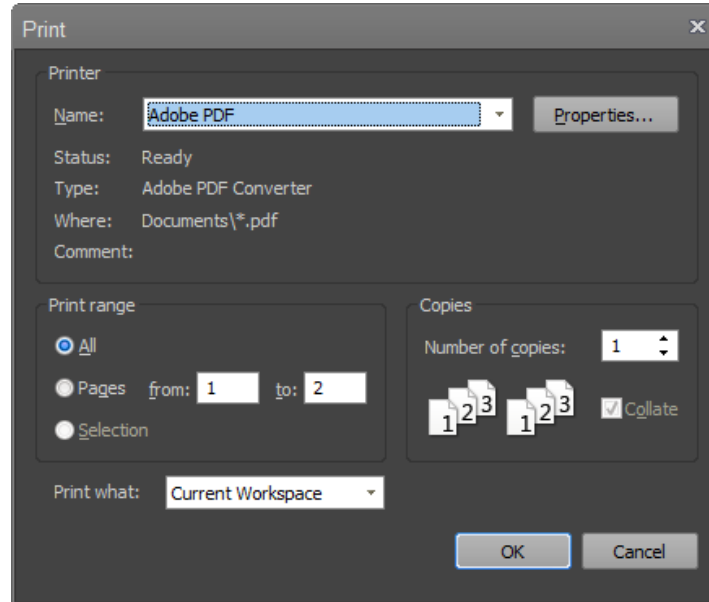
3. To save the changes and close the dialog, click **OK**.  
To close the dialog without saving the changes, click **Cancel** or **X**.

## Print dialog

The **Print dialog** is a modified **Windows™ Print dialog**. All the controls on this dialog, except the **Print what** and **Print range**, are standard Windows™ functions and are not described here.

### Open the dialog

On the **File Ribbon tab** (“File tab” on page 72), select **File ▶ Print** or press **Ctrl+P** on the keyboard.



### Print range

This is the standard print range control. However, the **Print range** options are disabled when the **Print dialog** is opened from the **Experiment Explorer context menu ▶ Batch Print dialog** (“Batch Print dialog” on page 734). In this case, **All** option is selected and the entire print queue is processed.



## Print what options

- The **Print what** options are not displayed when the **Print dialog** is opened from the **Experiment Explorer context menu** ▶ **Batch Print dialog** (“Batch Print dialog” on page 734).
- If the **Print dialog** is opened from any of the standard application views (Workspace, Heat Map, Results, or Overlays), the **Print what dropdown** displays the following options:
  - **Current Workspace**: Prints the currently active **Workspace** in the order specified in the **Page setup dialog** (“Page Setup dialog” on page 728).
  - **Heat Map**: Prints the **Heat Map** and its legend scaled to one page width.
  - **Results**: Prints **Results tables** (both primary and secondary, if available) scaled to one page width.
  - **Overlays**: Prints **Overlay plots** and their corresponding legends (if enabled) with one plot per row. The **Overlay galleries** cannot be printed.
  - **Sample List**: Prints the **Sample List** as described in “Print the Sample List” (“Sample List context menus” on page 227).
- If the **Print dialog** is opened from the **Performance Test module** (Chapter 20, “Performance Test reports”), the **Print what dropdown** list displays the following options:
  - **Selected PT Report**: Prints the currently active and visible **Performance Test report** (“Current PT results” on page 561). If no PT report is visible, the most recent PT report is printed.
  - **Levey-Jennings**: Prints the currently selected **Levey-Jennings Report** for all installed detectors. Only the selected date range and statistics are printed. If no Levey-Jennings is selected, the most recent Levey-Jennings plot is printed using the specified date criteria.
  - **Performance History**: Prints the **performance history** for the currently selected baseline. If no baseline is selected, the most recent performance history is printed.
- Selecting one of the options changes the view to the selected view and sends the selected option to the print queue when the **OK** button is clicked.
- If any selection is not present or does not contain data, nothing is printed.
- The default selection for the **Print what dropdown** is determined by the view currently in focus.

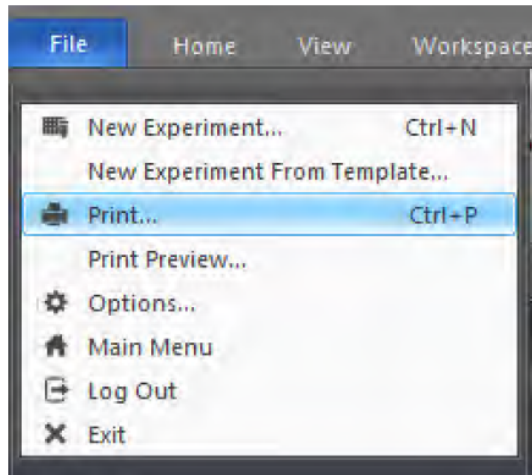
## Print Preview dialog

The **Print Preview dialog** displays the document as it will appear when printed and lets you set various printing options.

### Open the dialog

On the **File Ribbon tab** (“File tab” on page 72), select **File ▶ Print Preview**. The **Print Preview** option is enabled only when the instrument is not acquiring.

The **Print Preview dialog** opens as a full screen modal dialog that displays the document as it will appear when printed.



## Print Preview dialog

The controls to close the print preview, print the document, select portrait vs. landscape, open the page setup dialog, view the document full page width, view a whole page or multiple pages, and navigate are displayed in a command bar at the top of the dialog.

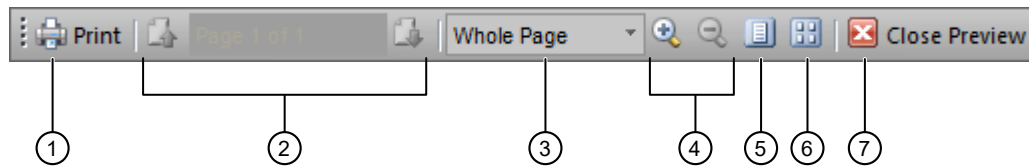


Figure 139 Print Preview dialog controls

- |   |                        |
|---|------------------------|
| ① Print   | ④ Zoom In and Zoom Out |
| ② Page Up and Page Down controls and Current page | ⑤ Page Size            |
| ③ Page View dropdown                              | ⑥ View Multiple Pages  |
|   | ⑦ Close Preview        |

- To print the document, click **Print**, which opens a **Print dialog** that contains standard Windows™ print functions.
- To navigate through the preview pages, click the **Page Up** and **Page Down** controls. The current page number is displayed in the **Current page** box. Alternatively, navigate directly to a specific page by entering the page number in the **Current page** box, then press the **Enter** key.
- To resize the page view, you can select any of the following options:
  - Select a size from the **Page View** dropdown. By default, whole page is selected.
  - **Zoom In** and **Zoom Out** controls allow the page view to zoom in and out. When the page zoom is at 500%, **Zoom In** is disabled. When the page zoom is at 10%, **Zoom Out** is disabled.
  - Click the **Page Size** button to set the view to the whole page.
  - To view up to 9 pages at one time, click the **View Multiple Pages** button, then select the pages to view from the menu.

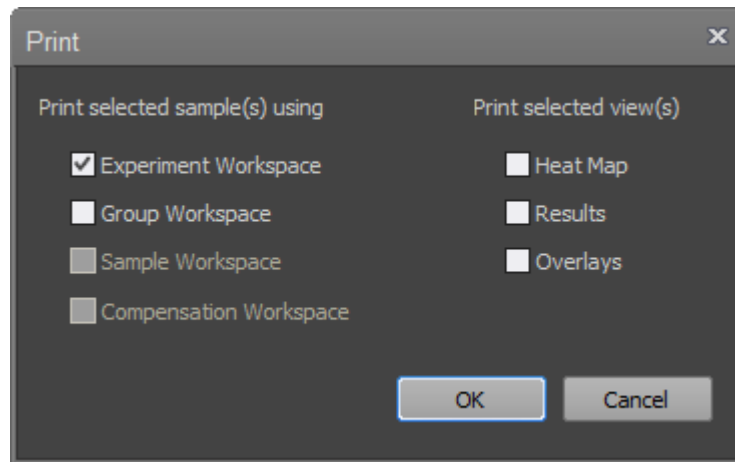


- To close the **Print Preview dialog**, click **Close Preview**.

## Batch Print dialog

### Open the dialog

The **Print dialog** is displayed when the **Print** option is selected from the **Experiment**, **Group**, or **Sample context menu** in the Experiment Explorer or from the **Heat Map context menu**.



### Print dialog options

#### Print selected sample(s) using

- The options in this group allow you to select the Workspaces to use when printing the selected samples. Available options are:
  - **Experiment Workspace**
  - **Group Workspace**
  - **Sample Workspace**
- You can select multiple Workspaces. By default, the **Experiment Workspace** is selected.

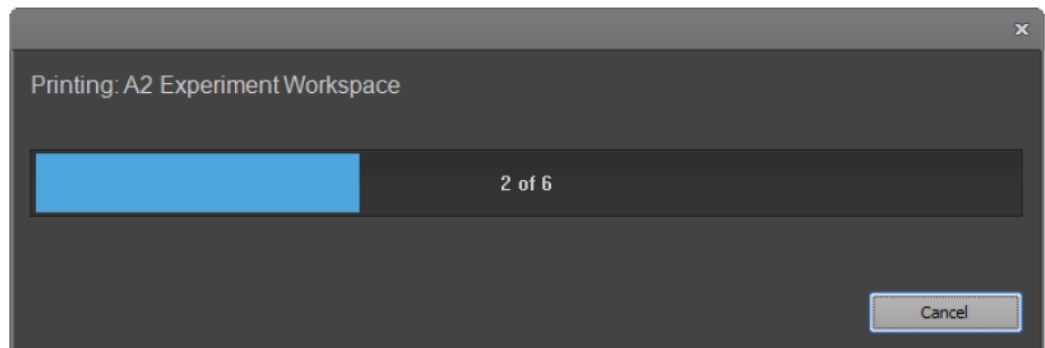
#### Print selected view(s)

- The options in this group allow you to select the view to print. Available options are:
  - **Heat Map**
  - **Overlay**
  - **Results**

By default, none of these options are selected.

- To start printing, click **OK**, which opens the **Print dialog** (“Print dialog” on page 730).
- Before printing, the **Heat Map**, **Overlays**, and **Results view** are refreshed to ensure that all statistics are up to date.

- During the batch print process, the progress of the print job is indicated on a dialog. The number of items in the print queue is the sum of Workspaces printed for each Sample and views selected to print.



- To cancel the print job, click **Cancel** on the progress dialog.

## Print order

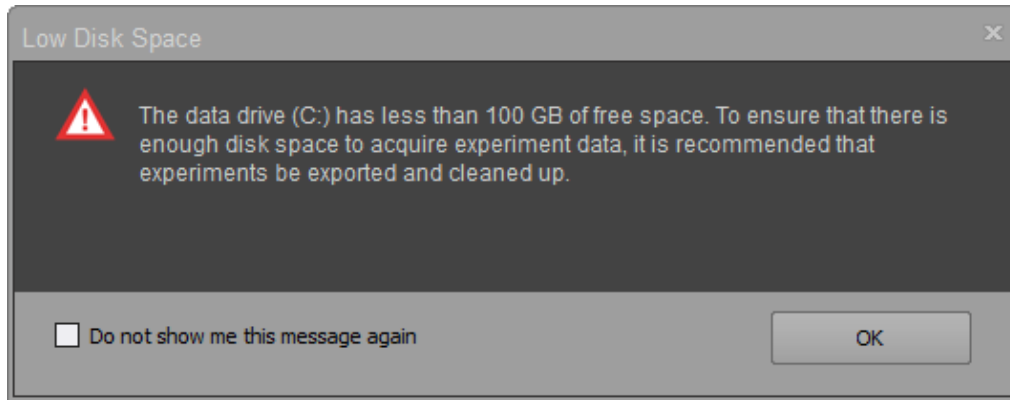
- The print order is:
  - Experiment Workspace (<n>)
  - Group Workspace (<n>)
  - Sample Workspace (<n>)
  - Results
  - Heat Map
  - Overlays

where <n> is the sample index

- A page break is added between each new Workspace and each new Sample, and between the Heat Map, Results, and Overlays.
- Items are printed using the current page layout settings.
- The Heat Maps legend is scaled to the available print width.
- Result tables are scaled to the page width.
- The Results view is based on the currently selected filters and statistics.
- Overlays are printed in the order that they appear on the Overlay view (Chapter 9, “Overlays”). Gallery plots are not printed.

## Low Disk Space dialog

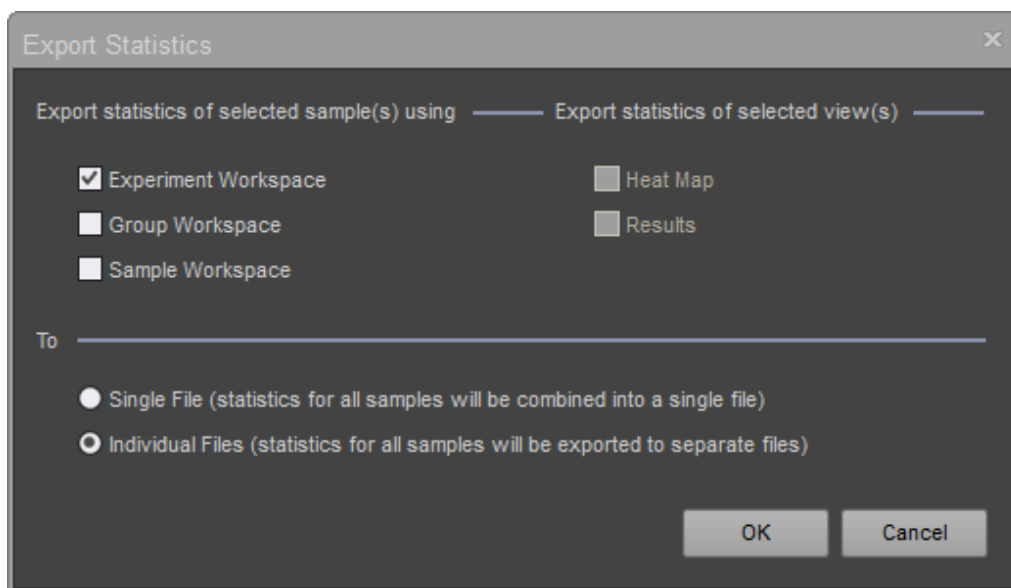
The **Low Disk Space** dialog is displayed when activating Experiments, if the **Warn on Low Disk Space** option is enabled in the **Administrator** tab of the **Options** dialog and the disk space is less than the specified threshold on the primary data (image data and experiment data) drives.



- By default, when activating experiments, the Attune™ Cytometric Software displays the **Low Disk Space** dialog if the available disk space is less than 100 GB on the primary data (image data and experiment data) drives.
- You can ignore the **Low Disk Space** warning and suppress the dialog for the length of the active session by selecting the **Do not show me this message again** option.
- You can disable the warning or change the low disk threshold in the **Administrator** tab of the **Options** dialog (see “Administrator” on page 676).

## Export Statistics dialog

The **Export Statistics** dialog is displayed when **Export Statistics** is selected from the **Plate**, **Sample**, **Group**, or **Experiment context menus** of the **Experiment Explorer** (“Experiment Explorer context menus” on page 301).



### Export statistics of selected samples using

The options in this group let you select the Workspaces to use when exporting the selected Samples. You can select multiple Workspaces. Available options are:

- **Experiment Workspace**
- **Group Workspace**
- **Sample Workspace**

By default, the **Experiment Workspace** is selected. For Workspace statistics, the individual plot statistics are not exported.

### Exports statistics of selected views

The options in this group let you export the statistics of selected views. You can select multiple views. Available options are:

- **Heat Map**
- **Results**

By default, the no views are selected.

## To

The options in this group allow you to select the output. Available options are:

- **Single File:** Exports the statistics as a single CSV file. Results are collated into a single table for the Workspace statistics and the **Results** tab's primary table.
- **Individual Files:** Exports the statistics as multiple CSV files. The software creates a CSV file for each workspace and each selected sample. If the **Heat Map** and **Results** views are selected, the software creates extra files.
- To start exporting, click **OK**. When prompted, complete the:
  - **Save** dialog (“File Save (Export) dialog” on page 715) for the **Single File** option.
  - **Open** dialog (“File Open (Import) dialog” on page 721) for the **Individual Files** option.
- To close the dialog without printing, click **Cancel** or **X**.



## Derived gate dialog

The *Derived Gate dialog* allows you to create gates based on Boolean operators (i.e., *Derived gates*).

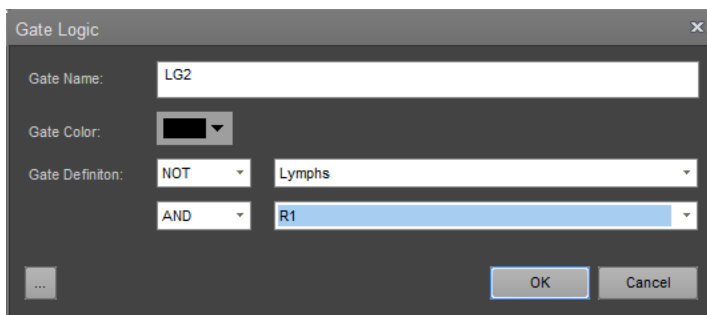
### Launch the dialog

In the Workspace tab (“Gating Tools group” on page 83), click **Derived Gate** to open the Derived Gate dialog.



The Derived Gate button is enabled when the Workspace view is displayed and the current Workspace has at least one gate present.

### Collapsed view

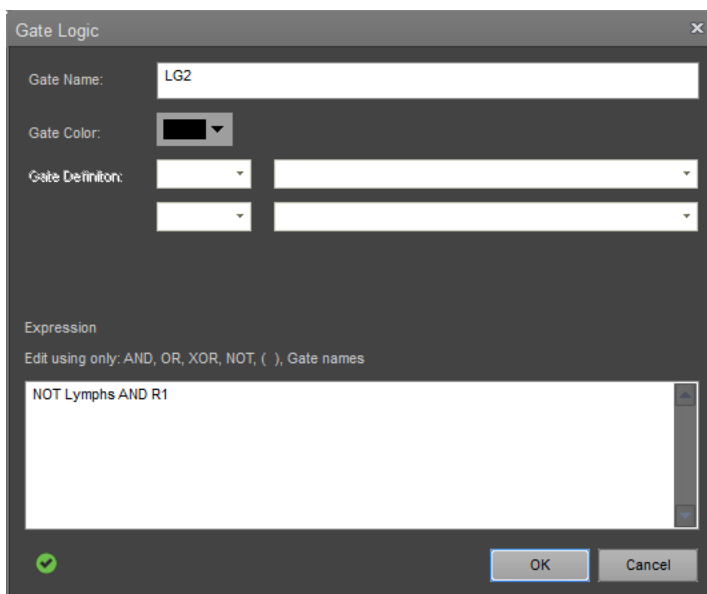


The Gate Logic dialog in collapsed view shows the following fields:

- Gate Name: LG2
- Gate Color: [Black]
- Gate Definiton: NOT (operator), Lymphs (gate)
- AND (operator), R1 (gate)

Buttons: OK, Cancel

### Expanded view



The Gate Logic dialog in expanded view shows the following fields:

- Gate Name: LG2
- Gate Color: [Black]
- Gate Definiton: [Empty]
- Expression: NOT Lymphs AND R1

Buttons: OK, Cancel

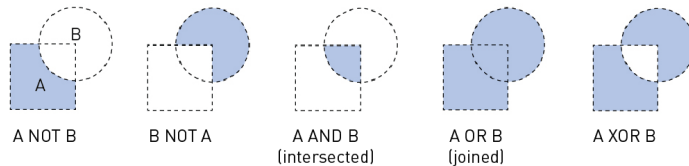
---

**Note:** The contents of the Derived Gate dialog changes depending on the context of the active Workspace.

---

## Boolean operators

The Boolean operators available for creating Derived gates are blank, AND, OR, XOR, AND NOT, and OR NOT.



## Derived gate dialog

### Gate name

Allows you to accept the default gate name, or enter a new name.

Gate Name:

- You can enter up to 50 characters.
- Duplicate names and names that are only a whole number are not permitted.
- The default name for a gate is **LG1**. If a gate using the default name already exists, the next available integer is automatically used; for example, **LG2**.

### Gate color

Allows you to select a gate color from the dropdown menu.

Gate Color:

### Gate definition

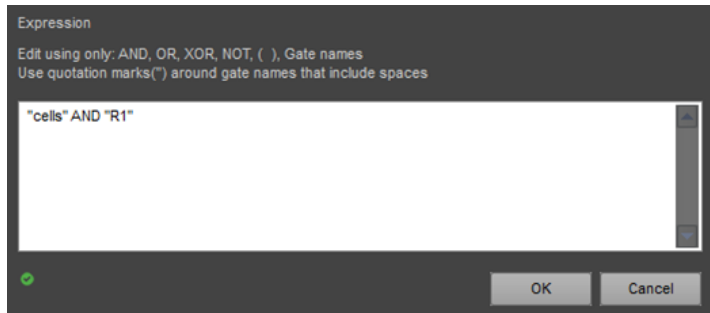
Allows you to create gate math expressions using Boolean operators.

Gate Definition:

- Select a gate from each of the two dropdown menus. The menus contain alphabetized lists of all gates on the active Workspace.
- From the first Boolean dropdown menu, select **<blank>** or **NOT**.
- From the second Boolean dropdown menu, select another Boolean operator **<blank>**, **AND**, **OR**, **XOR**, **AND NOT**, or **NOT**.

## Expression editor

- Click the **expander button** to expand the *Derived Gate dialog* and display the *Expression Editor*. If you made selections in the Gate Definition dropdown menus (above), they appear in the text box.



- Edit the text using only: **AND, OR, XOR, NOT, <blank>**, and gate names. You can enter at least 1500 characters.
- All gate names must be entered in quotation marks (i.e., enter R1 as “R1”).
- As you enter text, the software validates the expression and indicates the status:
  - ✓ = valid
  - ⚠ = invalid
- To save the changes and close the dialog, click **OK**.
- To close the dialog without creating gates, click **Cancel** or **X**.

## Edit gates dialog

The *Edit Gates* dialog provides a list of all available gates in the active Workspace, and allows you to:

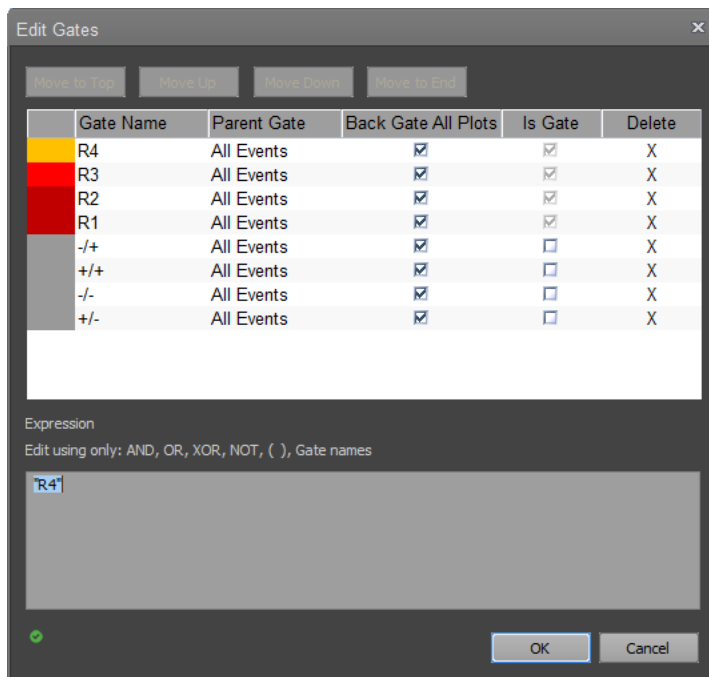
- Edit the gate color, gate math expression, and the order in which the gates are displayed in the dialog
- Delete gates

### Launch the dialog

In the Workspace tab (“View tab” on page 77), click **Edit Gates**.



The Edit Gates button is available when the Workspace view is displayed and the current Workspace has at least one gate present.



---

**Note:** The contents of the *Edit Gates* dialog changes, depending on the context of the active workspace.

---

## Edit gates dialog

### General properties

- By default, gates are added to the top of the list.
- Changes are applied to all gates after **OK** is clicked.
- To remove the gate name from the list and delete the gate from the plot, click **Delete**.
- You cannot sort the table or rearrange the columns.
- You can resize columns by selecting and dragging the column divider or by double clicking on the divider to auto-fit the column width to the contents.
- Vertical and horizontal scrollbars are displayed as necessary.

### Reorder gates

- To reorder the gate list, select a gate, then drag and drop it to a different location within the list. Changing the gate list order does not change the gating hierarchy.
- Alternatively, select a gate, then click **Move to Top**, **Move Up**, **Move Down**, or **Move to End**, as necessary.



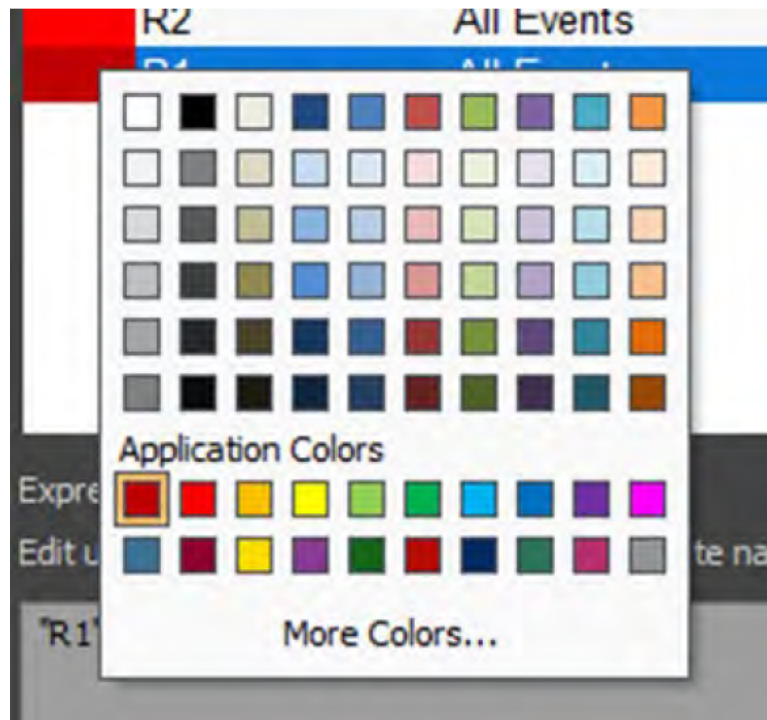
- To select multiple gates at one time, press **Ctrl** or **Shift** when selecting the gates. The order of the selected gates is maintained when they are moved.

## Resize columns

- To resize a column, drag the column divider, or double-click the divider to auto-fit the column width to the contents.

## Change gate color

- To change the gate color, select the **Color** cell to open the **color picker**.



## Edit gate names

- To edit a gate name, select a name from the dropdown menu in the **Gate Name** cell, or double-click the gate name and enter new text.
- You can enter up to 50 characters.
- Duplicate names and names that are only a whole number are not permitted.
- The default name for a gate is LG1. If a gate using the default name already exists, the next available integer is automatically used; for example, LG2.

## Backgate all plots

- The **Backgate all plots** option is checked by default for all newly created oval, rectangle, and polygon gates, if the **Backgate all plots** option is checked in the Gate options dialog ("Gate Options" on page 656).
- When selected, backgating is applied in all dot and precedence density plots for the selected gate. This paints the events contained within the gate on these plot types based on the color of the gate.
- When de-selected, backgating is turned off for the selected gate on all plots.

## Is gate

- The “Is Gate” column as a checkbox control indicates whether a region has an associated gate.
- By default, histogram, oval, rectangle, and polygon regions are gates. Derived gates are always gates.
- When a region is associated with a gate, it cannot be removed and the checkbox is disabled.
- The software limits the number of total gates to 128. Because each quadrant contains four members, quadrants are created initially without an associated gate. For quadrant members that do not have an associated gate, the “Is Gate” checkbox is unchecked. When the “Is Gate” is checked for a quadrant member, a gate is created for that quadrant member and the checkbox is disabled.

## Edit gate equation

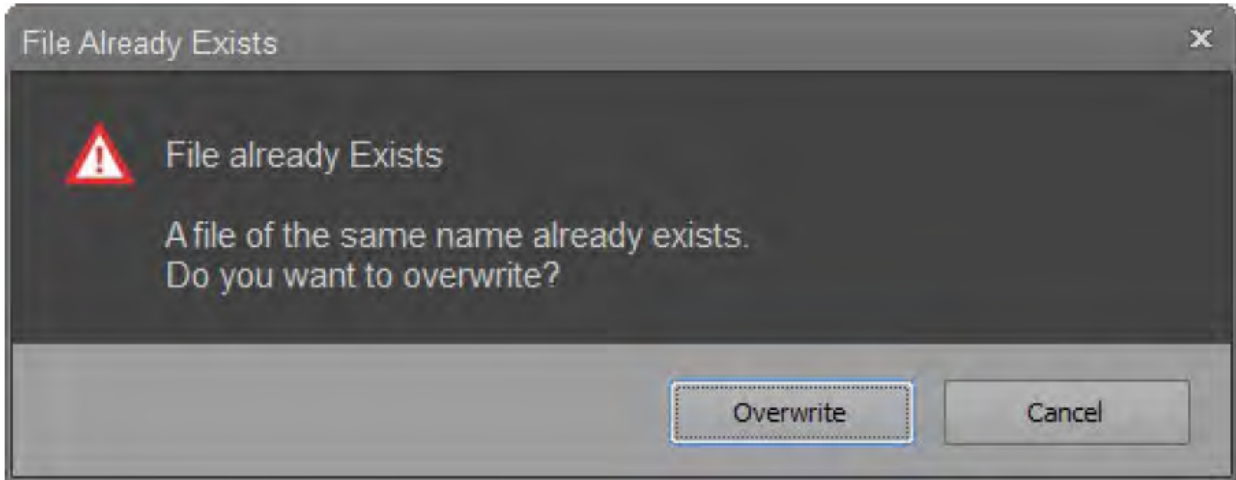
- To edit a gate equation, select a derived gate, then enter the equation in the *Expression Editor*.



- Only derived gate equations can be edited; otherwise this feature is disabled. If you select a non-derived gate, the Expression Editor is disabled.
- Edit the text using only: **AND, OR, XOR, NOT, <blank>**, and gate names. Gates names must be bounded by quotation marks (“gate name”). You can enter at least 1500 characters.
- As you enter text, the software validates the expression and indicates the status:
  - ✓ = valid
  - ⚠ = invalid
- To save the changes and close the dialog, click **OK**.

## Confirm file save dialog

The *Confirm File Save* dialog helps to prevent exported files from being accidentally overwritten. The dialog is automatically displayed when you attempt to export a file to a selected export location that already contains a file of the same name.

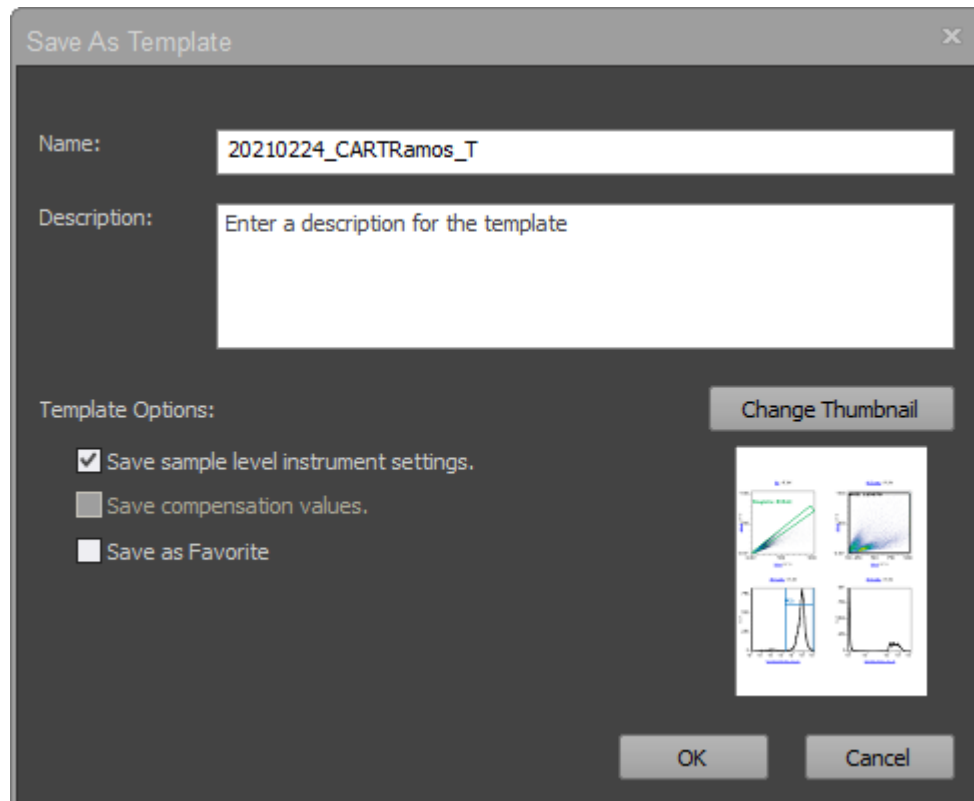


- To replace an existing file with the current file, click **Overwrite**.
- If you are exporting a batch of files, and multiple duplicate file names exist in the export location, the **Do this for all other cases (x found)** check box is enabled, where **x** is the number of existing files.
- To overwrite all duplicate files in the export location, check the **Do this for all other cases (x found)** check box.
- To decide on a case-by-case basis for each file, leave the check box unchecked.
- The checked state applies only to the current export batch and is not persisted.
- To close the dialog without exporting any files, click **Cancel** or **X**.



## Save As Template dialog

The **Save As Template** dialog enables you to save the current Experiment as a template so that the Experiment Workspace, Experiment Instrument Settings, Run Protocols, Heat Map Settings, and Experiment hierarchy are preserved for future Experiments.



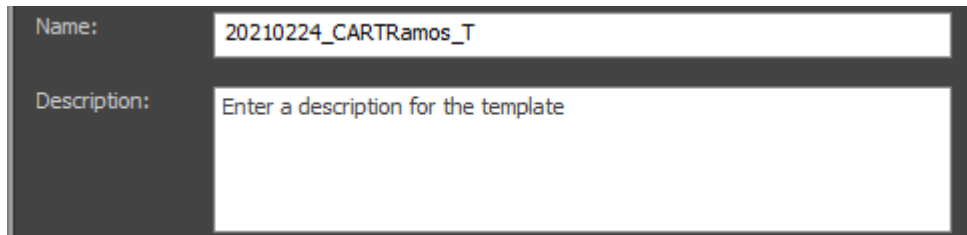
### Open the Save As Template dialog

Click the **Save As Template** icon in the **Home** tab (“Home tab” on page 74) or use the keyboard shortcut **Ctrl+S**. This saves the current experiment as a template.



Alternatively, right-click a **Tube** or **Plate Experiment**, then select the **Save As Template** option.

## Name and Description fields

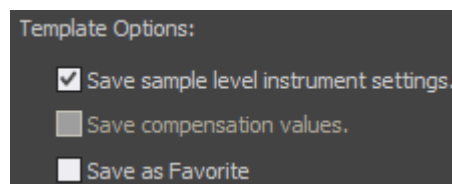


The screenshot shows a dialog box with two input fields. The first field is labeled "Name:" and contains the text "20210224\_CARTRamos\_T". The second field is labeled "Description:" and contains the placeholder text "Enter a description for the template".

- By default, the **Name** text field shows the name of the current Experiment from which the **Save As Template** is selected, with the name of the Experiment appended with “\_T”. If the default name is already in use, the name is checked for the suffix “(#)”, where # is the next available integer. The template name field allows up to 50 printable characters.
- The **Description** field is optional and allows up to 500 printable characters.
- If the template name is in use when **OK** is pressed, a warning dialog indicates that the template already exists.
  - **Overwrite** saves the template and replaces the existing template.
  - **Cancel** closes the dialog and returns to the **Save as Template** dialog.

## Template options

There are three options to save an Experiment as a Template:

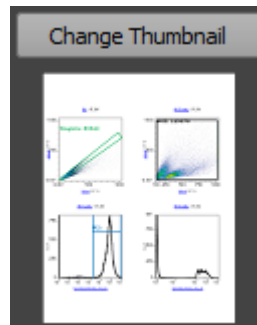


The screenshot shows a dialog box titled "Template Options:" with three checkboxes. The first checkbox, "Save sample level instrument settings.", is checked. The second checkbox, "Save compensation values.", is unchecked. The third checkbox, "Save as Favorite", is unchecked.

- **Save sample level instrument settings** is the default option and specifies whether the template is saved with the sample level instrument settings.
- **Save compensation values** option enables you to save the template with the compensation settings values (compensation XML). The option is deselected by default. If the experiment does not have compensation settings, this option is inactive. If a plate experiment contains Compensation controls mapped to wells, these wells are removed when this option is selected. If the option is unchecked, the template maintains the compensation controls to be used in the template, but the compensation settings values are not carried over.
- **Save as Favorite** specifies that the template is set as a favorite. This option is unchecked by default. Templates that are marked as “Favorite” are tagged for easy identification in the Momentum™ Software when using the Attune™ NxT instrument integrated with the Orbitor™ RS2 or the Orbitor™ RS3 Microplate Mover.

## Template thumbnail

Displays a thumbnail image that corresponds to the selected template.



- By default, the template thumbnail is the first page of the Experiment Workspace of the Experiment.
- You can select another template thumbnail by clicking **Change Thumbnail**, which opens the **File Open** browser and enables you to select images of the following formats: JPG, GIF, BMP, PNG, TIF, EMF.

## Template attributes

A template is saved with the following attributes:

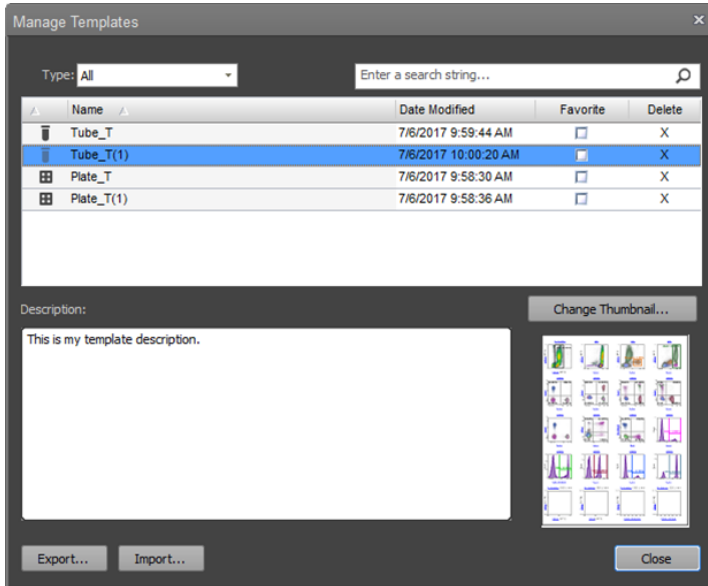
- Experiment hierarchy/layout (numbers of Groups and Samples, types of Samples)
- Heat Map mapping (location of Tubes and/or Wells)
- Workspaces (Experiment, Group and Sample)
- Instrument settings (Sample-level are optional, if selected as described above)
- Run Protocols
- Compensation settings (optional, as described above)
- Heat Map settings
- Overlay settings
- Results View settings (selected statistics, sorting, grouping, column sizing, column order)
- Experiment keywords
- Sample level keyword values

## Manage templates dialog

*Manage Templates* dialog allows you to edit, import, export, and delete templates. It contains the following elements:

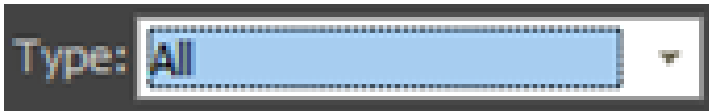
- Type filter (“Type filter” on page 750)
- Keyword search (“Keyword search” on page 750)
- Template list (“Template list” on page 751)
- Description field (“Description” on page 752)

- Thumbnails and Change Thumbnail button (“Thumbnails” on page 752)
- Export button (“Export” on page 754)
- Import button (“Import” on page 754)



## Type filter

Allows you to filter the templates by **type**. By default, the type filter is set to “All”.



- When **All** is selected, all templates that are valid for the active instrument model (i.e., BRVY) are displayed.
- When the filter type is changed to **Tube**, only tube templates are displayed.
- When the filter type is changed to **Plate**, only plate templates are displayed.

## Keyword search

Allows you to apply a search keyword to filter the list of displayed templates, where the experiment name or description contains the keyword.



- To filter the list of displayed templates, enter a keyword, then press **Enter** or click the **Search** button. This applies the search keyword as a filter, so that only the templates that contain the keyword in the name or the description field are displayed in the list.
- When a search keyword is applied, the search button displays “Clear”.
- To remove the keyword filter, click **Clear**, or delete the text in the edit control and press **Enter**.

## Template list

The template list is populated with the list of templates based on the active model and the filters applied by the experiment type and keyword search filters.

Name	Date Modified	Favorite	Delete
Tube_T	7/6/2017 9:59:44 AM	<input type="checkbox"/>	X
Tube_T(1)	7/6/2017 10:00:20 AM	<input checked="" type="checkbox"/>	X
Plate_T	7/6/2017 9:58:30 AM	<input type="checkbox"/>	X
Plate_T(1)	7/6/2017 9:58:36 AM	<input type="checkbox"/>	X

- The template list has the following columns:
  - **Template type:** Displays an icon that corresponds to the experiment type (tube icon for tube experiment and plate icon for plate experiment).
  - **Template name:** Displays the name of the template.  
To edit a template name, double-click on the template name. Alternatively, press the F2 key, if only a single template is selected.  
You can enter up to 50 alpha-numeric characters for the template name.  
The template name cannot end with a period character and cannot be any of the following words: “CON”, “PRN”, “AUX”, “CLOCK\$”, “NUL”, “COM1”, “COM2”, “COM3”, “COM4”, “COM5”, “COM6”, “COM7”, “COM8”, “COM9”, “LPT1”, “LPT2”, “LPT3”, “LPT4”, “LPT5”, “LPT6”, “LPT7”, “LPT8”, “LPT9”.
  - **Date Modified:** Displays the date the template was created or last modified. The date is displayed in system locale in the long time format (i.e., 6/15/2009 1:45:30 PM)
  - **Favorite:** Favorite checkboxes allow you to designate a template as a favorite.
  - **Delete:** Deletes the template.
- To sort a column, click on the desired column heading. When sorting, a sort arrow indicates the sort direction (i.e., ascending or descending). Favorite and Delete columns are not sortable.
- To select a template, click on the row that displays the template you wish to select. You can select multiple templates.
- When a single template is selected in the template list, the Description field updates to display the template description (if entered) and the Thumbnail field displays the corresponding thumbnail. If multiple templates are selected, the description field and thumbnails are not displayed.
- Right-click a template to open the template list context menu with options to **Export** or **Delete** the template.

## Description

Displays the optional, user-defined template description (see “Save As Template dialog” on page 747).

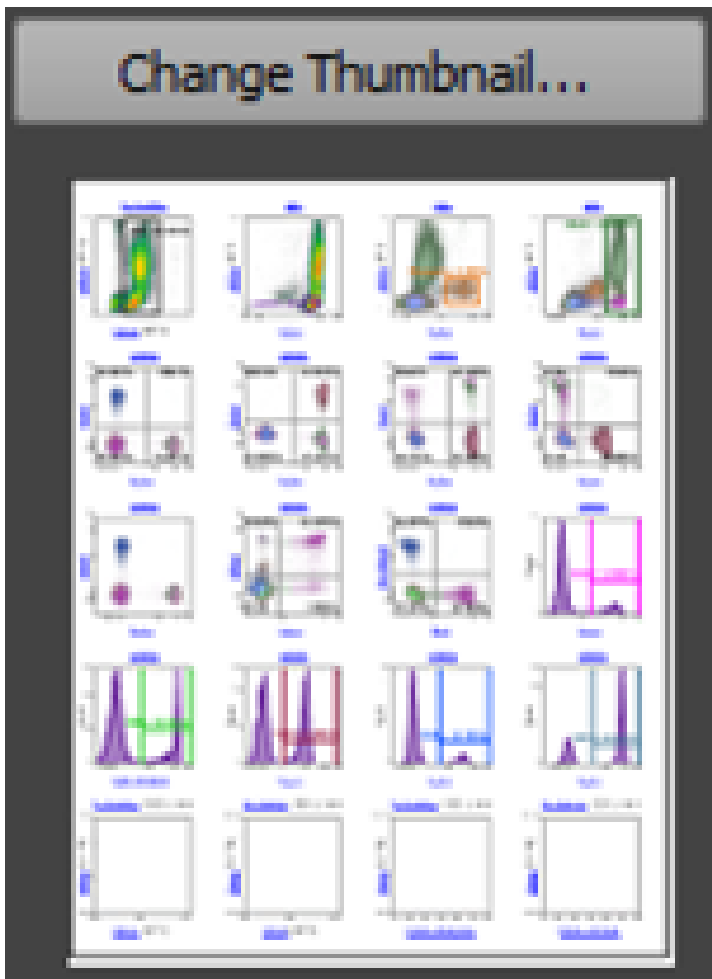


To edit the template description, click on the Description field and type in the desired description. You can enter up to 500 alpha-numeric characters.

## Thumbnails

Displays a thumbnail image that corresponds to the selected template.

- By default, the template thumbnail is the first page of the experiment workspace of the experiment.
- You can select another template thumbnail by clicking the **Change Thumbnail** button, which opens the File Open browser and allows you to select images of the following formats: JPG, GIF, BMP, PNG, TIF, EMF.



## Export

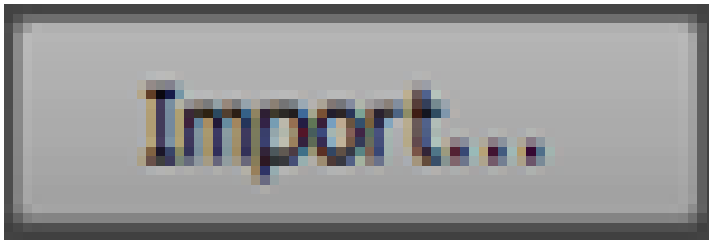
Allows you to export the selected dialog with your changes.



- Clicking **Export** when only a single template is selected opens the *File Save (Export) dialog* (“File Save (Export) dialog” on page 715) with the “File name” set to the selected template name and the “Save as type” set to Attune™ Experiment Template (\*.aet).  
Click **Save** on the File Save (Export) dialog to save the selected template to the specified location.
- Clicking **Export** when multiple templates are selected opens the *Folder Browser dialog* (“Folder Browser dialog” on page 723).  
Click **Select Folder** on the Folder Browser dialog to exports the selected templates to the selected folder with each template named as it was named in the Template list and given the .aet extension.

## Import

Allows you to import an Attune™ Experiment Template (\*.aet) file to your list of templates.



- Clicking **Import** opens the *File Open (Import) dialog* (“File Open (Import) dialog” on page 721) with the filter set to Attune™ Experiment Template (\*.aet).
- Select a \*.aet file to add the template to your list of templates and select the template in the Template list.
- You can import multiple templates. If you import multiple templates, the last imported template is selected in the Template list.



## Delete

- Click the **X** in the Template list or right-click a template and select **Delete** from the context menu to delete the selected template.

When you click **X** or select the context menu **Delete** option, the software displays the *Confirm Delete dialog*.

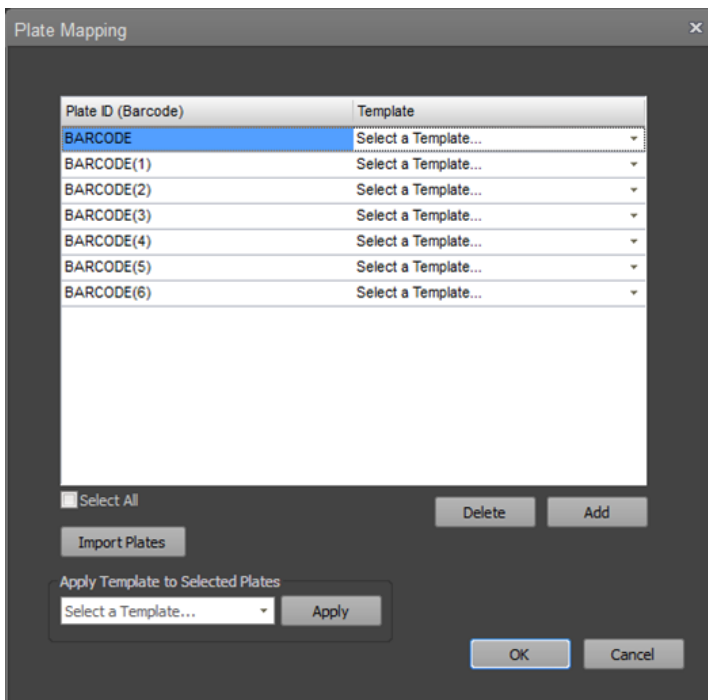
The dialog contains an option for **Do this for all other selected items (n found)**, where n is the number of remaining templates.

- Click **Yes** to delete the first selected template. The dialog will be presented again for each remaining selected template.
- Check the **Do this for all other selected items (n found)**, then click **Yes** to delete all selected templates.
- Click **Cancel** to return to the Manage Templates dialog without deleting any of the selected templates.

## Plate Mapping dialog

The **Plate Mapping** dialog enables you to map templates to specific plates that are defined by barcodes in the **Automation mode**. The **Plate ID (Barcode)** to **Template** mappings are accessible by the automation software (such as the Thermo Scientific™ Momentum™ Software) to run the selection of plates.

To open the **Plate Mapping** dialog, click the **Map Plates** button in the **Instrument ribbon tab** (“Automation group” on page 94).



The Plate Mapping dialog contains the following controls:

- **Plate ID (Barcode) to Template table** (“Plate ID (Barcode) to Template table” on page 757)
- **Select All** checkbox (“Select All” on page 757)
- **Delete** and **Add** buttons (“Delete and Add buttons” on page 758)
- **Import Plates** button (“Import Plates” on page 758)
- **Apply Template to Selected Plates** (“Apply Template to Selected Plates” on page 758)
- **OK** and **Cancel** buttons (“OK and Cancel buttons” on page 759)

## Plate ID (Barcode) to Template table

The **Plate ID (Barcode) to Template** table enables you to select a **template** from a dropdown list to assign to the selected **Plate ID** (i.e., barcode).

Plate ID (Barcode)	Template
BARCODE	Select a Template...
BARCODE(1)	Select a Template...
BARCODE(2)	Select a Template...
BARCODE(3)	Select a Template...
BARCODE(4)	Select a Template...
BARCODE(5)	Select a Template...
BARCODE(6)	Select a Template...

Select All Delete Add

- The **Plate ID (Barcode) to Template** table has these columns:
  - **Plate ID (Barcode):** Displays the Plate ID (Barcode). By default, the Plate ID name is **BARCODE**.  
If the name is already in use, the name is checked for the suffix “(#)”, where # is the next available integer. For example, BARCODE(2), BARCODE(3) etc.  
To edit a **Plate ID** name, double-click on the name. Alternatively, press the **F2** key, if only a single template is selected.  
You can enter up to 50 alpha-numeric characters for the template name.  
The template name cannot end with a period character and cannot be any of the following words: “CON”, “PRN”, “AUX”, “CLOCK\$”, “NUL”, “COM1”, “COM2”, “COM3”, “COM4”, “COM5”, “COM6”, “COM7”, “COM8”, “COM9”, “LPT1”, “LPT2”, “LPT3”, “LPT4”, “LPT5”, “LPT6”, “LPT7”, “LPT8”, “LPT9”.
  - **Template:** Enables you to select a **template** to map to a **Plate ID (Barcode)**.  
To map a **template** to a selected **Plate ID**, click **Select a Template...**, then select from the list of templates that have been marked as favorites.  
Alternatively, click **Select a Template...**, then select **Import/Browse...** at the top of the dropdown list, which opens the **Select Template** dialog (“Select template dialog” on page 760).

## Select All

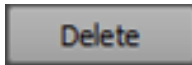
**Select All** selects all **Plate ID/Template** rows in the **Plate ID (Barcode) to Template** table.



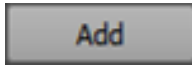
Deselect the checkbox to deselect all rows except the last row in table.

## Delete and Add buttons

- **Delete:** Deletes all selected rows from the **Plate ID (Barcode) to Template** table.

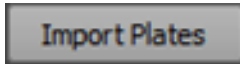


- **Add:** Adds a new row to the **Plate ID (Barcode) to Template** table using the default Plate ID name.

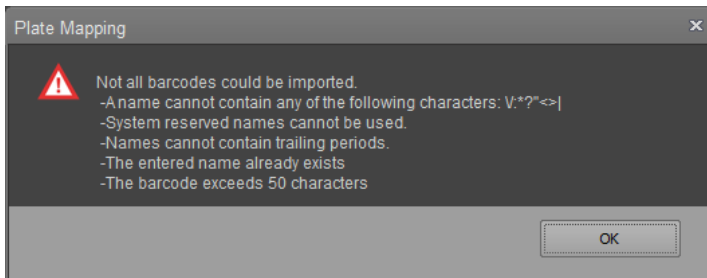


## Import Plates

**Import Plates** opens the **File Open (Import)** dialog (“File Open (Import) dialog” on page 721), which enables you to import a CSV file to add new barcodes to the **Plate ID (Barcode) to Template** table.

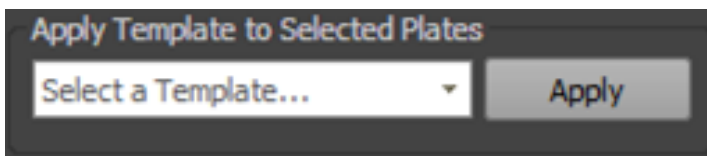


The CSV file is read from column by column and row by row and only names that are permitted are added. If an invalid name is detected, a dialog is displayed stating the reason that not all barcode names were imported:



## Apply Template to Selected Plates

**Apply Template to Selected Plates** enables you to assign a template to the selected **Plate ID/Template** rows.



- Click **Select a Template...** to select from the list of templates that have been marked as favorites. Alternatively, click **Select a Template...**, then select **Import/Browse...** at the top of the dropdown list, which opens the **Select Template** dialog (“Select template dialog” on page 760).
- Click **Apply** to apply the selected template to the selected rows.

## OK and Cancel buttons

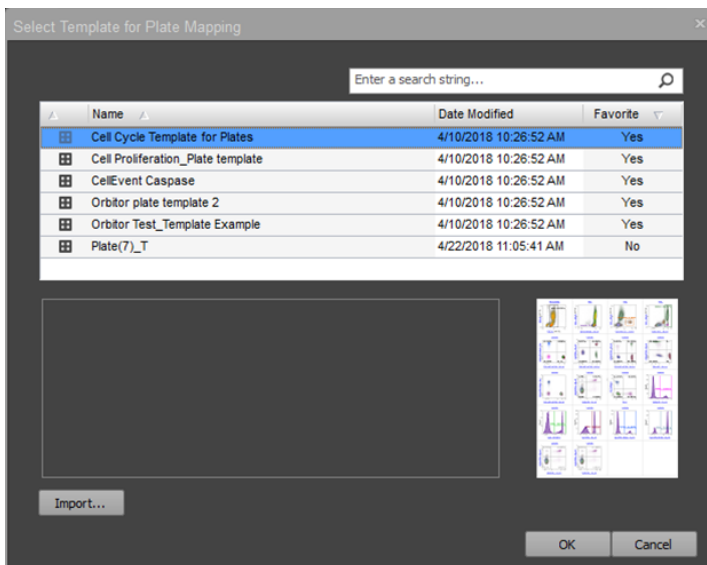
- Click **OK** to save the **Plate ID (Barcode) to Template** mappings in the database. The **Plate ID (Barcode) to Template** mappings are accessible by the automation software (such as the Thermo Scientific™ Momentum™ Software) to run the selection of plates.
- Click **Cancel** to close the dialog without saving the mappings.

## Select template dialog

The *Select Template* dialog is a modal dialog that allows you to select a Template to map to the selected Plate IDs (Barcodes).

The Select Template dialog only shows Plate templates for the active instrument model (e.g., BRVY). Tube templates are NOT displayed.

- Template Keyword search (“Select template dialog” on page 760)
- Template list (“Template list” on page 761)
- Template description (“Template description” on page 761)
- Thumbnails (“Thumbnails” on page 762)
- Import (“Import” on page 762)
- OK and Cancel (“OK and cancel buttons” on page 763)



### Template keyword search




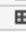
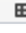

Allows you to apply a search keyword to filter the list of templates displayed in the template list, where the experiment name or description contains the keyword.



- To filter the list of displayed templates, enter a keyword, then press **Enter** or click the **Search** button. This applies the search keyword as a filter, so that only the templates that contain the keyword in the name or the description field are displayed in the list.
- When a search keyword is applied, the search button displays “X”.
- To remove the keyword filter, press the **X** button, or delete the text in the edit control and press **Enter**.

## Template list

The Template list is populated with a list of templates based on the active model and the filters applied by the Experiment type and keyword search filters.

Name	Date Modified	Favorite
 Cell Cycle Template for Plates	4/10/2018 10:26:52 AM	Yes
 Cell Proliferation_Plate template	4/10/2018 10:26:52 AM	Yes
 CellEvent Caspase	4/10/2018 10:26:52 AM	Yes
 Orbitor plate template 2	4/10/2018 10:26:52 AM	Yes
 Orbitor Test_Template Example	4/10/2018 10:26:52 AM	Yes
 Plate(7)_T	4/22/2018 11:05:41 AM	No

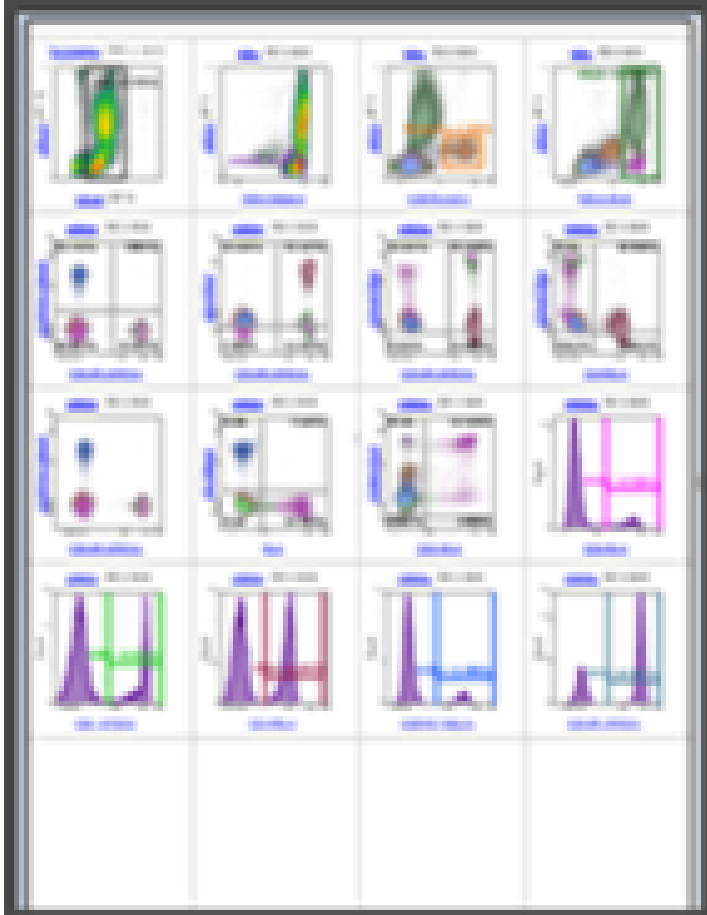
- The Template list has the following columns:
  - **Template type:** Displays an icon that corresponds to the Experiment type (tube icon for Tube experiment and plate icon for Plate experiment). Only Plate experiment templates are listed in the “Select Template for Plate Mapping” dialogue.
  - **Template name:** Displays the name of the template.
  - **Date modified:** Displays the date the template was created or last modified. The date is displayed in system locale in the long time format (i.e., 6/15/2009 1:45:30 PM)
  - **Favorite:** Shows whether the template was designated as a favorite.
- To sort a column, click on the desired column heading. When sorting, a sort arrow indicates the sort direction (i.e., ascending or descending). Favorite and Delete columns are not sortable.
- To select a template, click on the row for the template you wish to select. You can select multiple templates.
- When a single template is selected in the template list, the Description field updates to display the template description (if entered) and the Thumbnail field displays the corresponding thumbnail. If multiple templates are selected, the description field and thumbnails are not displayed.
- Right-click a template to open the template list context menu with options to **Export** or **Delete** the template.

## Template description

Displays the optional, user-defined template description (see “Save As Template dialog” on page 747).

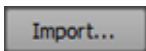
## Thumbnails

Displays a thumbnail image that corresponds to the selected template. By default, the template thumbnail is the first page of the experiment workspace of the experiment.



## Import

Enables you to import an Attune™ Experiment Template (\*.aet) file to your list of templates.



- Click **Import** to open the **File Open (Import)** dialog (“File Open (Import) dialog” on page 721) with the filter set to Attune™ Experiment Template (\*.aet).
- Select a **\*.aet file** to add the template to your list of templates and select the template in the **Template list**.
- You can import multiple templates. If you import multiple templates, the last imported template is selected in the **Template list**.



## OK and cancel buttons

- Click **OK** to close the dialog and assign the selected template to the Template list that has launched the dialog.
- Click **Cancel** to close the dialog. The template list in the Plate Mapping dialog reverts to its previously mapped template. The default is blank.

## Map Sample List Data dialog

The **Map Sample List Data** dialog enables you to create Samples and map them based on location when you import a **Sample List** from a CSV file.

**Note:** \*.xls files can not be imported as **Sample Lists**.

Sample Information	Column
Location (\$WELLID)	LOCATION
SAMPLE (\$FIL)	SAMPLE
SAMPLE_NOTES (\$CELLS)	SAMPLE_NOTES
GROUP (\$SMNO)	GROUP
GROUP_NOTES (\$SRC)	GROUP_NOTES
EXPERIMENT (\$PROJ)	EXPERIMENT
EXP_NOTES (\$COM)	EXP_NOTES
PLATE (\$PLATENAME)	PLATE
PLATE_ID (\$PLATEID)	PLATE_ID
PLATE_NOTES	PLATE_NOTES

Has header row

OK Cancel

- When the **Sample List** is imported, the software looks for the following default system-defined columns/keywords:
  - LOCATION
  - SAMPLE
  - SAMPLE\_NOTES
  - GROUP
  - GROUP\_NOTES
  - EXPERIMENT
  - EXP\_NOTES
  - PLATE
  - PLATE\_ID
  - PLATE\_NOTES
- The **Map Sample List** dialog also displays the keyword label and the corresponding control for each custom (user-defined) keyword.

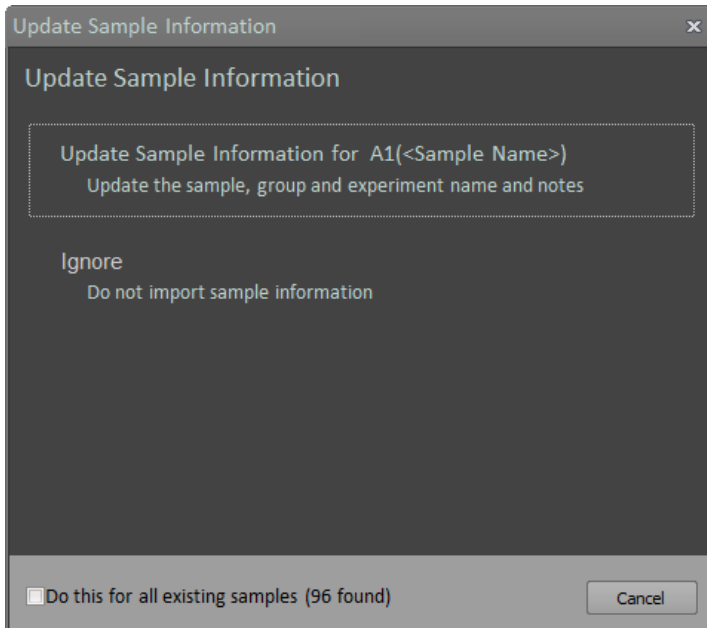
- The added keywords mapped from a CSV file are based on the existing Experiment keywords. If there are no custom keywords assigned in the Experiment, the dialog only displays the system-defined keywords and default sample list columns, regardless of the custom keywords present in the CSV file.
- If the column headers match the specified column names, they are mapped to their respective fields.
- The **Has header row** option enables the first row in the CSV file to be interpreted as a column header instead of actual row data. By default, this option is selected.



- When checked, the first row of the CSV is interpreted as column names instead of actual row data, and the combo box controls in the dialog are populated with the column header values as parsed from the CSV file.
  - When unchecked, the first row is interpreted as actual row data, and the combo box controls are populated with generic column names (Column 1, Column2, etc.) with the number of columns corresponding to the number of columns in the CSV file.
  - The combo box controls also include a **None** option, which excludes the corresponding sample information from being read from the CSV file.
- Click **OK** to update the Experiment with the Sample List data in the CSV file using the column mapping specified in the dialog.  
Click **Cancel** or **X** to close the dialog without updating the Experiment and importing the Sample List data.

## Update sample information dialog

The **Update Sample Information** dialog is displayed if you attempt to import a Sample List from an existing CSV file, but Samples already exist in the Experiment or Plate (i.e., Samples are assigned to a location).

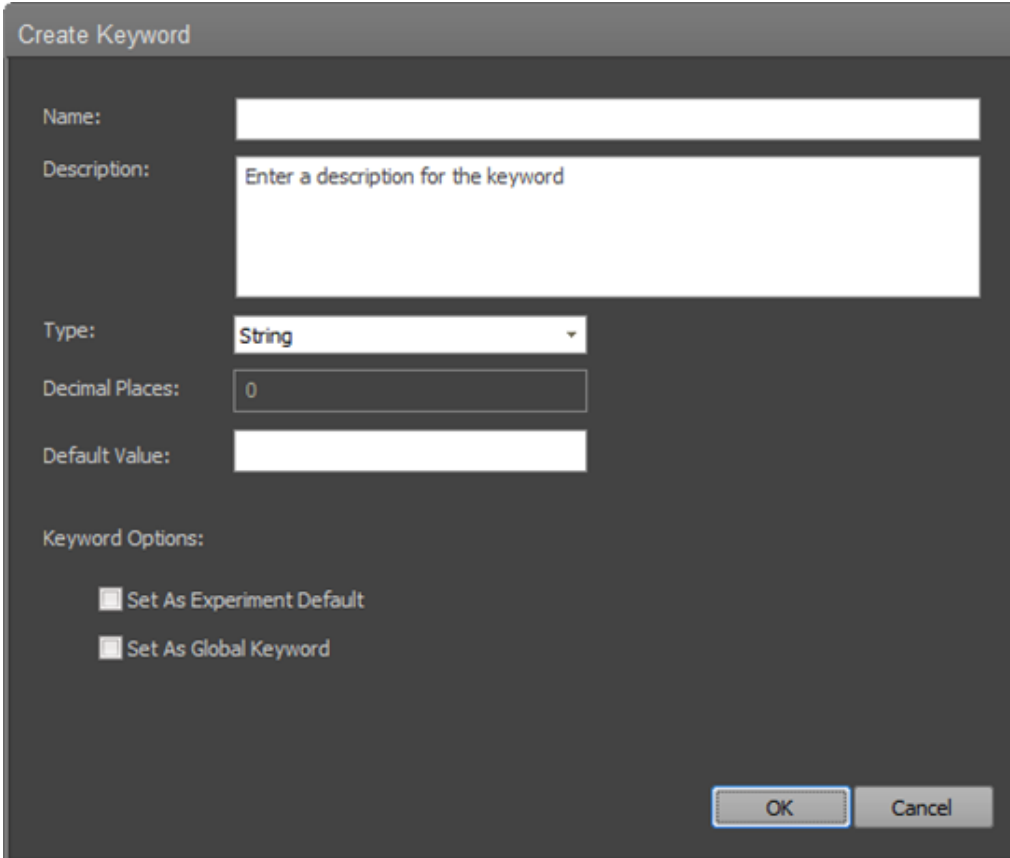


- Click **Update Sample Information** to update the Sample, Sample Notes, Group, Group Notes, Experiment, Experiment Notes, Plate, Plate ID, and Plate Notes for the specified Sample.
- Click **Ignore** to skip the import/update of the Sample information for the specified Sample.
- Check **Do this for all existing samples** to repeat the Update or Ignore action for every Sample in the Experiment or Plate.
- Click **Cancel** or the **X** button to close the dialog and prevent additional actions from being performed (any changes to existing Samples before this are not reverted).

## Create/Edit keyword dialog

The **Create/Edit Keyword** dialog enables you to define new keywords to add to the global keyword options list (see “Plate Options” on page 675).

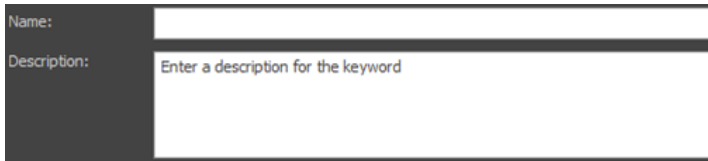
- When the dialog is launched from **Keyword options** by clicking the **Create** button, the dialog title is **Create Keyword**.
- When the dialog is launched from **Keyword options** by clicking the **Edit** button, the dialog title is **Edit Keyword**.



The screenshot shows a dialog box titled "Create Keyword" with a dark gray background. It contains the following fields and options:

- Name:** A text input field.
- Description:** A text area with the placeholder text "Enter a description for the keyword".
- Type:** A dropdown menu currently set to "String".
- Decimal Places:** A text input field containing the value "0".
- Default Value:** A text input field.
- Keyword Options:** Two checkboxes:
  - Set As Experiment Default
  - Set As Global Keyword
- Buttons:** "OK" and "Cancel" buttons at the bottom right.

## Name and Description



Name:

Description:

- The **Name** field is required and allows up to 50 printable characters.
  - If the **Name** field is left blank when **OK** is clicked, a warning dialog prompts you to enter a name for the keyword.
  - If the keyword name is in use when **OK** is clicked, a warning dialog indicates that the keyword name already exists.
- The **Description** field is optional and allows up to 500 printable characters.

## Type

**Type** dropdown enables you to select **String** or **Number** for the **keyword type**.

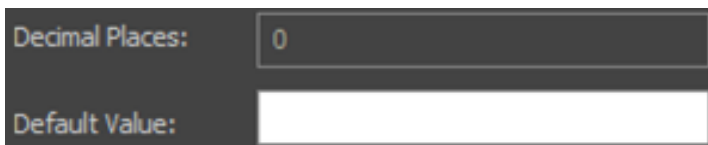


Type:

- If the **keyword type** is changed from **String** to **Number** or vice versa, a warning dialog prompts you to confirm the change.
  - Click **Yes** to confirm the change, close the warning dialog and the **Edit Keywords** dialog, and reset the keyword values.
  - Click **No** to close the warning dialog and return to **Edit Keywords** dialog.

## Default Value and Decimal places

**Default Value** enables you set the keyword value automatically when creating Samples in an Experiment using the selected keyword.



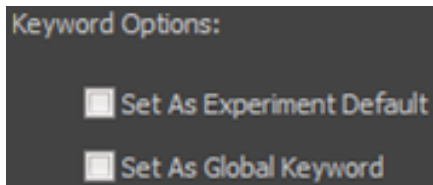
Decimal Places:

Default Value:

- The **Default Value** is blank by default and can be left blank.
- The **Default Value** field allows up to 50 characters.
- When the **Type** is set as a **Number**, the **Default Value** only accepts the negative sign, numbers, thousands separators, and decimal places as per the locale setting.
- If the **Type** is set to **Numeric**, the **Decimal Places** field is enabled. The **Decimal Places** setting is used to format a numeric keyword for display in the **Sample List** table and sets the number of decimal places to include in the FCS file.
  - The default number of decimal places is set to **0**.
  - The number of decimal places allows up to 30 decimals.

## Keyword Options

**Keyword Options** allow the keyword to be set as an **Experiment Default** and as **Global Keyword**.

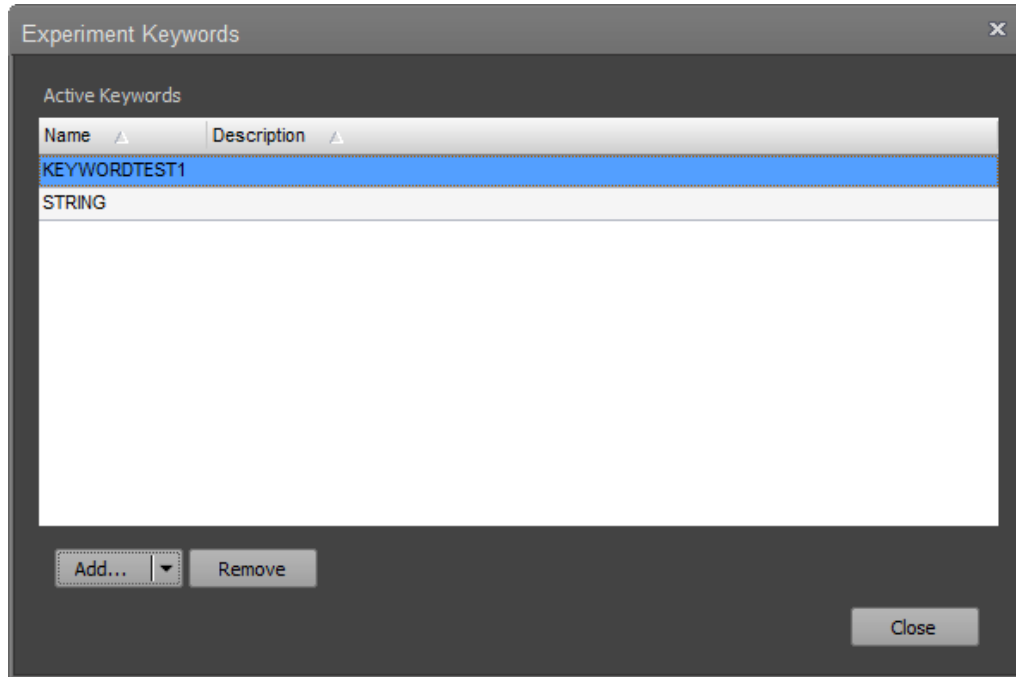


- **Set As Experiment Default** adds the selected keyword to the Experiment keyword list automatically when the Experiment is created.  
All users can set a keyword as an Experiment Default.
- **Set As Global Keyword** sets the custom keywords in the **Keyword Options** dialog list to **Global**.  
Global keywords are included in each user account after creation.  
The **Set As Global Keyword** checkbox is only visible to Administrator accounts, and only they can edit or delete a global keyword. User-level accounts cannot delete a global keyword.

## Experiment Keywords dialog

The **Experiment Keywords** dialog displays the **Active Keywords** list for the current Experiment and enables you to add, edit, and remove custom keywords for the Experiment.

The **Active Keywords** list has two columns: **Name** and **Description**.



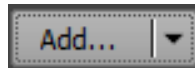
---

**Note:** The **Experiment Keywords** (including any local modifications made) persist in the Experiment. To add new keywords to the master list to be available for use in an Experiment, use the **Keyword tab** of the **Options** dialog (“Plate Options” on page 675) to edit or create user and global keywords.

---

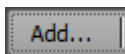
### Add/Edit

The **Add/Edit** split button enables you to add and edit custom keywords for the active Experiment.

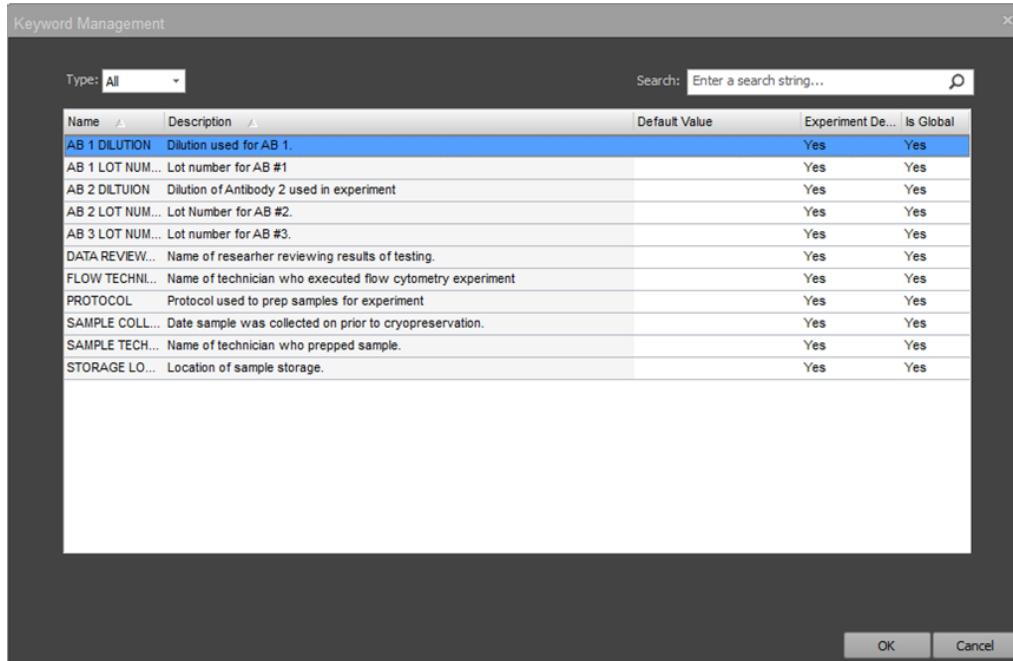


### Add

- Click **Add** to open the **Select Keywords** dialog, which includes all the keywords defined in the account that are not already included in the current Experiment.







- Keywords that are already part of the **Experiment Keyword** set are not displayed in the **Select Keywords** dialog.
- You can select multiple keywords, then click **OK** to copy them to the Experiment and show them in the **Experiment Keywords** dialog.
- Click **Cancel** or **X** to close the dialog without taking action.

## Edit

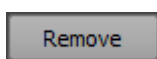
- Click the **Edit** part of the **Add/Edit** split button to open the **Edit Keyword** dialog described in “Create/Edit Keyword dialog” (“Create/Edit keyword dialog” on page 767).



- Changes made in the **Edit Keywords** dialog are only applied to the keywords in the current Experiment. Changes are not applied to the keyword of same name in the master list in the **Keyword tab** of the **Options** dialog (“Plate Options” on page 675).
- Click **OK** to apply the changes locally to the keyword in the current Experiment.
- Click **Cancel** or **X** to cancel any changes and close the dialog.

## Remove

- Click **Remove** to remove any selected keywords from the **Active Keywords** list of the current Experiment.

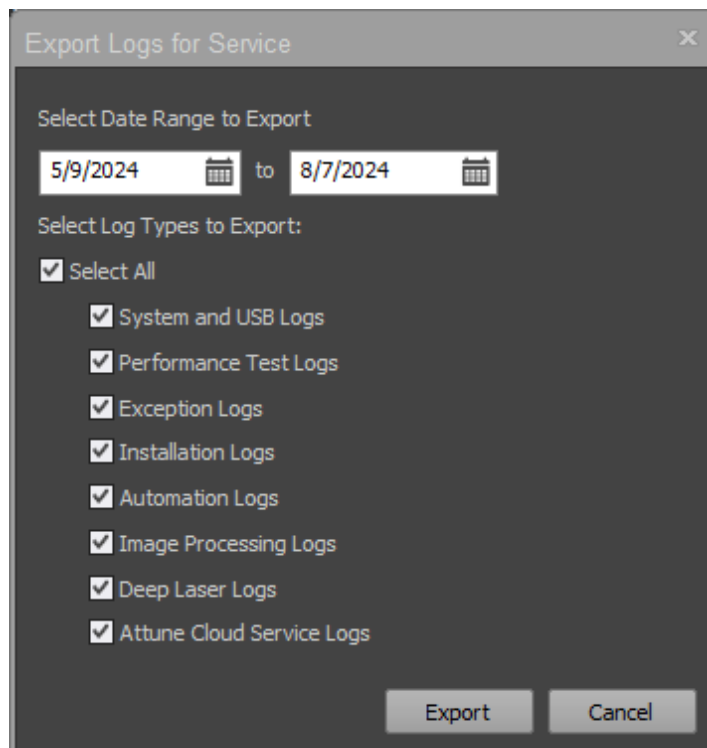


- You can select and remove multiple keywords together at once.
- Keywords that are removed from the **Active Keywords** list are not removed from the master list in the **Keyword tab** of the **Options** dialog (“Plate Options” on page 675).

## Export Logs for Service dialog

The **Export Logs for Service** dialog enables you to export selected logs to a single zip file based on a specified date range.

The dialog consists of two **date pickers** and a list of **log types** that you can select for export. By default, the **date range** is set to include the last 90 days of logs and all logs are selected.



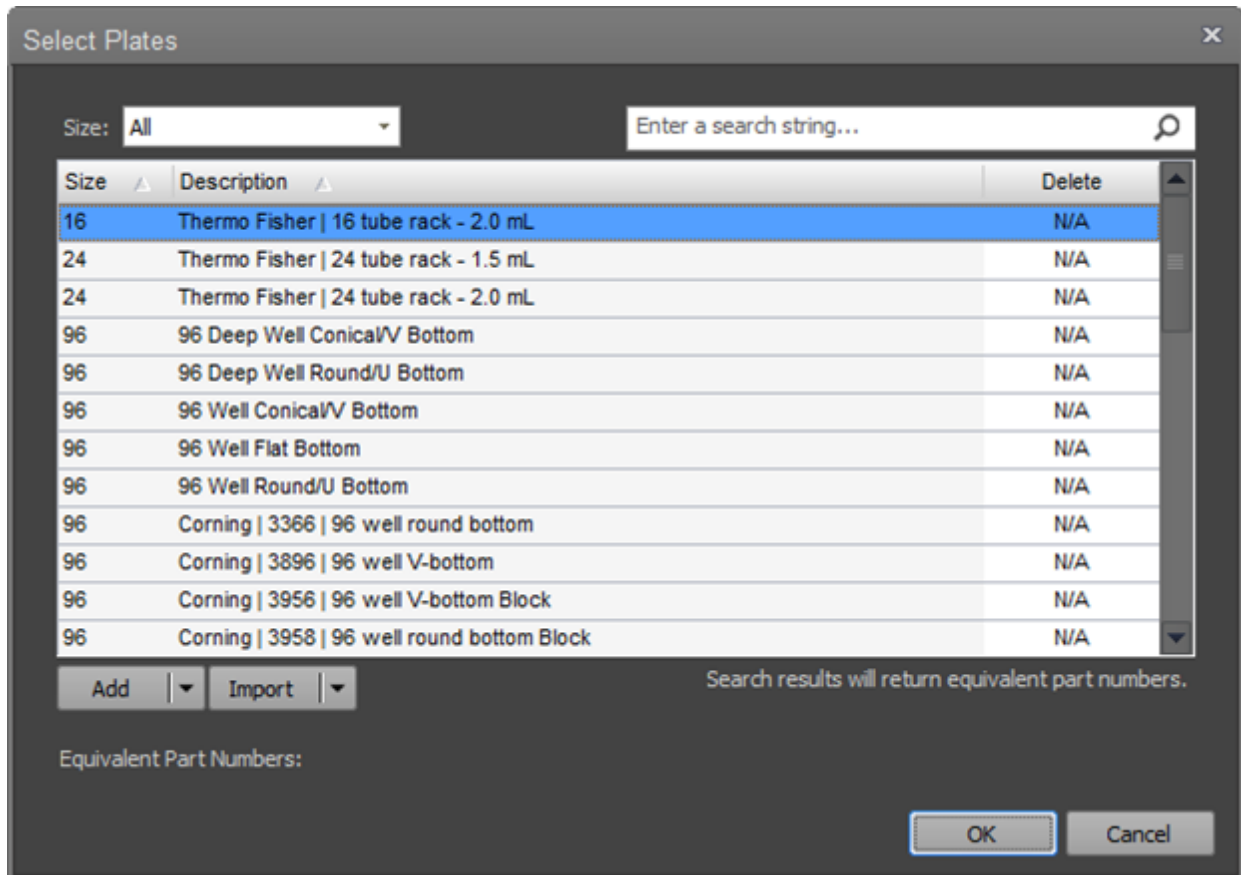
- **Select Date Range to Export** controls enables you to select the start and end dates. The start date must occur before the end date. If you enter an invalid date range, the instrument displays a warning.
- **Select Log Types to Export** enables you to export all logs at one time or select the individual log types you want to export.
  - **System and USB Logs** (which include the system and USB logs)
  - **Performance Test Logs** (which include the logs, corresponding iteration logs, and FCS files)
  - **Exception Logs** (only zipped logs are included)
  - **Installation Logs** (installation and firmware logs)
  - **Automation Logs**
  - **Image Processing Logs** (DMSService and other processing logs)
  - **Deep Laser Logs**
  - **Attune Cloud Service Logs** (logs for the service which connects the Attune™ Software to the Thermo Fisher Cloud)

- **Export** opens the **File Save (Export)** dialog (“File Save (Export) dialog” on page 715), which enables you to save the selected logs as a zip file in the desired folder.  
The default file name for the zip file is **AttuneServiceLogs\_YYYY-MM-DD\_HH-MM-SS**, where YYYYMMDD\_HHMMSS is the date and time.
- **Cancel** or the **X** button closes the dialog with no further action.

## Select Plates dialog

The **Select Plates** dialog displays a list of all available plates in the database. You can add these plates to the **Quick Select Plates** list in the **Options > Plate Options** dialog (“Plate Options” on page 675) or access them from the **New Experiment** dialog (“New Experiment dialog” on page 606) or the **Customize Plate Experiment** panel (“Customize experiment (Plate or Tube) options” on page 466).

**Note:** The **Select Plates** dialog is available only when a CytKick™ Max™ Autosampler is connected to the Attune™ instrument.

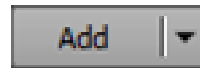


- The plates list has columns for **Size**, **Description**, and **Delete**. To sort the plates list by Size or Description, click the **Size** or **Description** column header.
- To search the available plates in the database, enter the search prompt in the **Search** field. You can search available plates based on **Size** or **Description**.
- When you perform a search, **equivalent part numbers** where the plate definitions agree with the original search prompt are displayed in the lower left of the dialog.

Equivalent Part Numbers:  
264574 | 264575 | 264576 | 264579 | 264675 | 264573

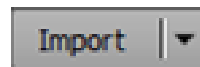
- To filter the available plate list by Size, select the desired option from the **Size** dropdown. Available options are **All**, **16**, **24**, **96**, and **384** (well plates).

- The **Delete** column is only visible for Administrator accounts when the dialog is launched from the **Quick Select Plates** list in the **Options ▶ Plate Options** dialog (“Plate Options” on page 675). When the mouse moves over the **Delete** column, the pointer changes to the hand pointer, if the plate can be deleted. You can only delete user-defined plates. System-defined plates show **N/A** in the **Delete** column and cannot be deleted. If a plate definition is in use as an existing Plate Experiment or Template, deletion is not allowed.
- To remove a Plate from the list of favorites, click the **X** in the **Delete** column next to the plate you want to delete.
- **Add/Edit/Duplicate** split button enables you to add, edit, or duplicate a plate definition.



Only Administrators, Advanced Users, and Service accounts can add, edit, or duplicate plate definitions. For users who do not have this permission, the button displays **View** and no dropdown list is provided.

- To add a new plate definition to the plates list, click **Add** to open the **Create Plate** dialog (“Create/Edit plate” on page 776).
- To edit a plate definition in plates list, select the **Edit** option on the **Add/Edit/Duplicate** split button to open the **Edit Plate** dialog (“Create/Edit plate” on page 776).
- To duplicate a plate definition, select the **Duplicate** option on the **Add/Edit/Duplicate** split button. The **Create Plate** dialog opens with all the fields prepopulated to agree with the selected fields in the plate definition. However, the description field is updated to have (#) appended to the description to create a unique name for the duplicated plate definition.
- **Import/Export** split button enables you to import or export a plate definition.



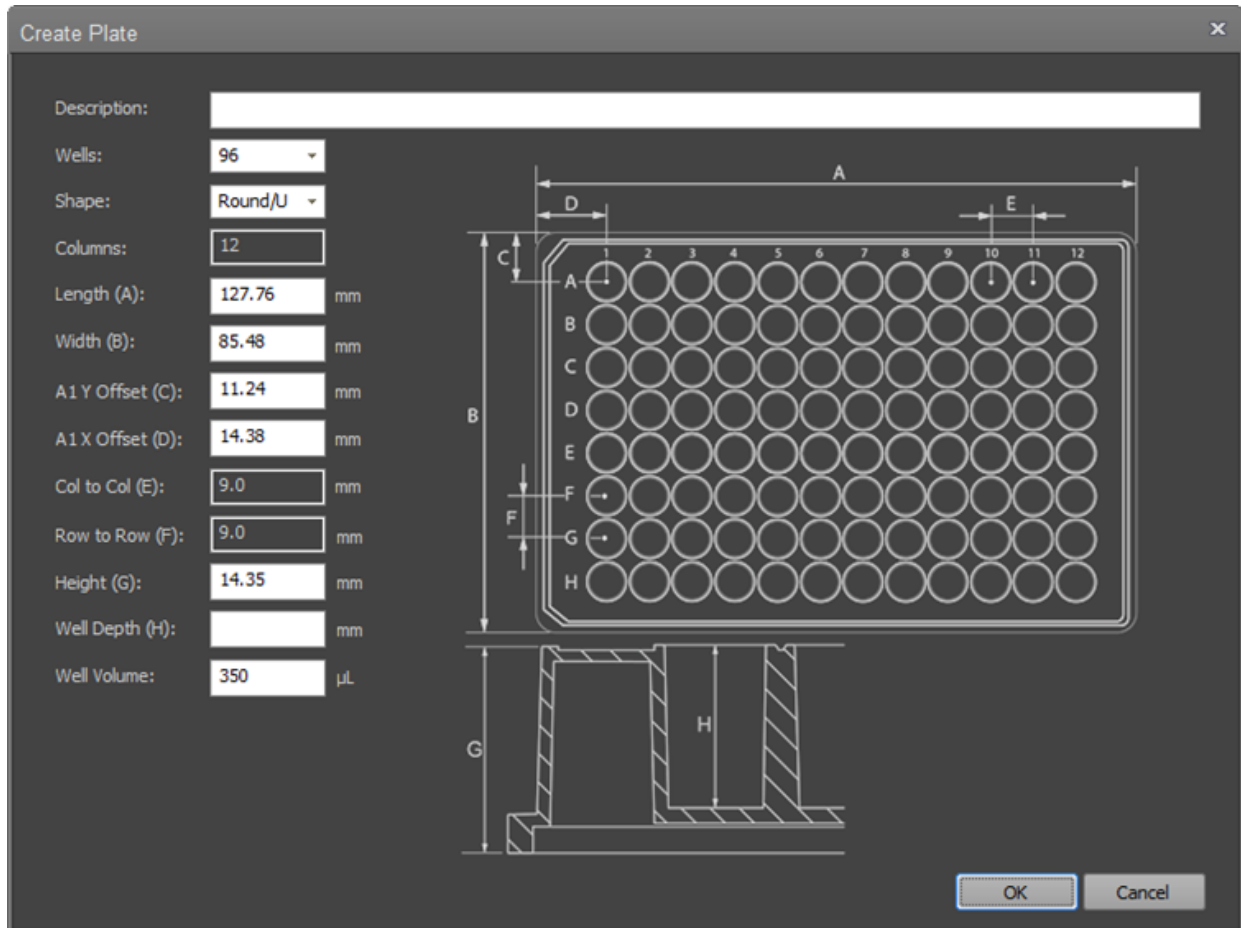
This button is a split button only when launched from the **Quick Select Plates** list in the **Options ▶ Plate Options** dialog (“Plate Options” on page 675); otherwise, it is a standard button with the **Import** option only.

- To import a plate definition to the plates list, click **Import**. The **File Open (Import)** dialog opens, which enables you to import a file containing plate definitions (“File Open (Import) dialog” on page 721).
- To export a plate definition, select the **Export** option on the **Import/Export** split button. The **File Save (Export)** dialog opens, which enables you to export the selected plate definitions (“File Save (Export) dialog” on page 715).

## Create/Edit plate

**Create/Edit Plate** dialog enables you to create custom plates with specific dimensions (such as plate height and well volume) to add to the database.

The **Create/Edit Plate** dialog is available only when a CytKick™ Max™ Autosampler is connected to the Attune™ instrument.



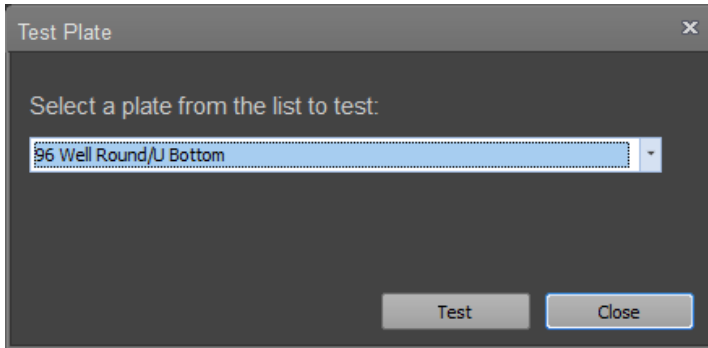
- If you have clicked **Add** in the **Select Plates** dialog (“Select Plates dialog” on page 774) to open the dialog, the dialog title is **Create Plate**.  
If you have clicked **Edit** in the **Select Plates** dialog to open the dialog, the dialog title is **Edit Plate**.
- If a plate is used in an existing Experiment or Template, the **Edit Plate** dialog does not allow change of any of the plate fields (they are read-only).
- The **Description** field must contain a unique description (it cannot be blank). The field allows 260 characters.
- New plate definitions have these default values:
  - **Wells:** 96 (available options are 48, 96, and 384 wells)
  - **Shape:** Round/U bottom (available options are Round/U, Conical/V, and Flat/F bottom)
  - **Columns (read-only):** 6 for 48-well, 12 for 96-well, 24 for 384-well
  - **Length (A):** 127.76 mm

- **Width (B):** 85.48 mm
- **A1 Y Offset (C):** 11.24 mm for 48-well and 96-well, 8.99 mm for 384-well
- **A1 X Offset (D):** 18.88 mm for 48-well, 14.38 mm for 96-well, 12.13 mm for 384-well
- **Col to Col (E):** 9 mm
- **Row to Row (F):** 9 mm
- **Height (G):** 14.35 mm
- **Well Depth (H):** blank
- **Well Volume:** 350  $\mu$ m

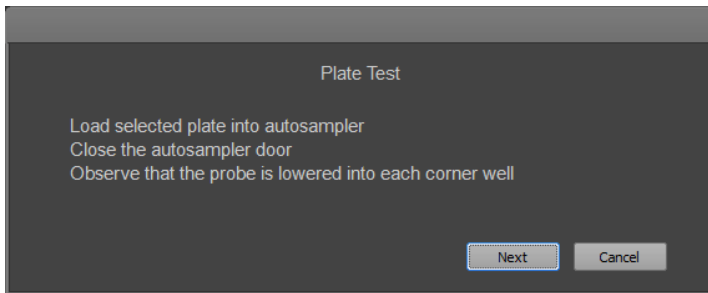
## Test Plate dialog

The **Test Plate** dialog enables you to validate a plate to ensure that the autosampler probe position is in the correct location in all four corners of the plate.

The dialog consists of a **Plate Selection** dropdown and **Test** and **Close** buttons.



- **Plate Selection** dropdown enables you to select a plate to test. The list of plates is based on the Quick Select plates definition.
- **Test** button initiates the test procedure and displays the test instructions.



## Run the plate test

1. Click **Test Plate** button on the **Instrument** ribbon tab ("Service group" on page 93) to open the **Test Plate** dialog.
2. Select the plate you want to test from the **Plate Selection** list, then click **Test**.
3. Load the selected plate into the autosampler, close the autosampler door, then click **Next**.



During the test, the selected plate definition is sent to the instrument. The autosampler then iterates through each corner of the selected plate or tube rack and lowers the probe to the well depth as defined in the plate definition. The autosampler cycles through each corner of the plate three times.

---

**Note:** You cannot cancel the Plate Test procedure after it has started.

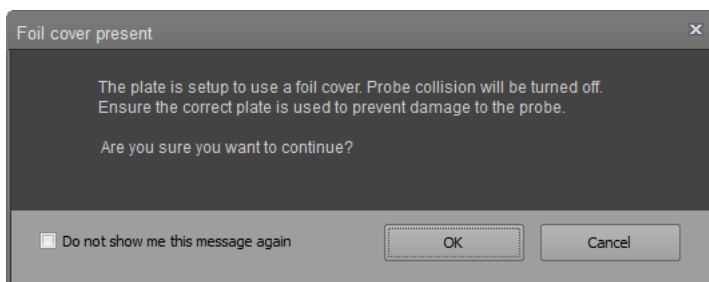
---



4. At the completion of the test, the dialog displays the message **Test completed without errors!**. If a probe collision error occurs, the error message is displayed instead of the successful completion message.
5. When the test is completed, click **Close** to close the **Test Plate** dialog.

## Foil cover present dialog

When **Use Foil Cover** is enabled (“Create a plate experiment” on page 609), the software displays the **Foil cover present** warning dialog at the beginning of a plate acquisition.

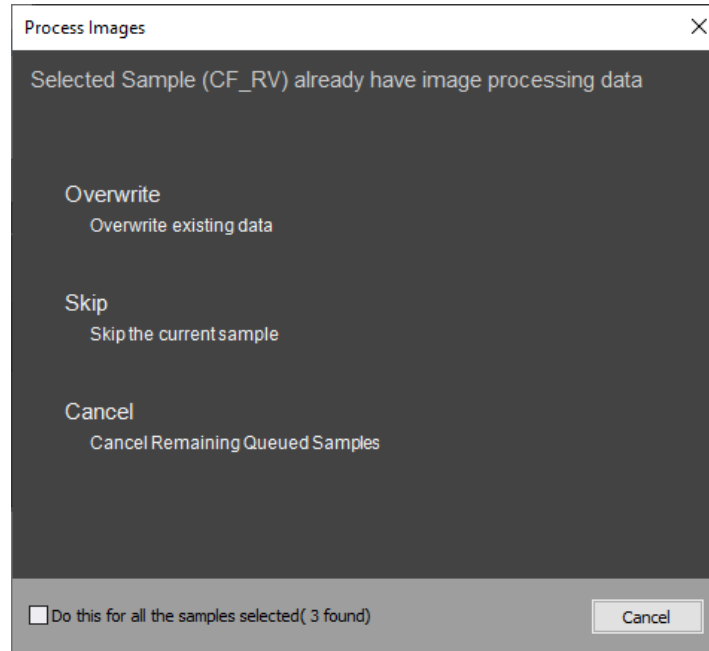


- The **Foil cover present** dialog is displayed when you click the **Setup Comp**, **Record Comp**, **Record All**, or **Record Plate** button on the **Collection Panel**, or if you attempt to **Run/Record** from a Manual well).  
The dialog is not displayed when you run a Tube sample or a Compensation sample from a Tube within a Plate Experiment.
- Click **OK** to close the dialog and proceed with acquisition.
- Click **Cancel** or the **X** button to abort the acquisition and close the dialog.
- Select **Do not show me this message again** to suppress dialog for the length of the current session.

## Process (Reprocess) Images dialog

The **Process (Reprocess) Images** dialog is displayed when you start image processing for a sample already has image processing data. The software displays the dialog for each sample in the processing request that meets this condition.

**Note:** You can reprocess images in a sample to allow a different processing model or algorithm, or to use other populations or features to extract cell image information. When samples are reprocessed, any ongoing processing is canceled and the new processing request is added to the queue.



The **Process (Reprocess) Images** dialog states that the **Selected Sample <sample name> already has image processing data**, and presents three options:

- **Overwrite:** Proceeds with the processing of the selected sample and overwrites the existing image processing data.
- **Skip:** Skips the samples with existing image processing data and processes only the samples that do not have existing data.
- **Cancel:** Cancels the processing request without adding the selected samples to the image processing queue. The image processing dialog remains open, allowing you to update the sample selection, if desired.

If you have selected multiple samples with existing image processing data, you can select **Do this for all the samples selected** to apply the action to all samples that meet this condition. When the last sample that meets this condition is confirmed, the image processing dialog is closed.

- If **Do this for all the samples** is selected and **Overwrite** is clicked, selected samples that already have image processing data are added to the queue and processed along with any samples that do not have existing data.

---

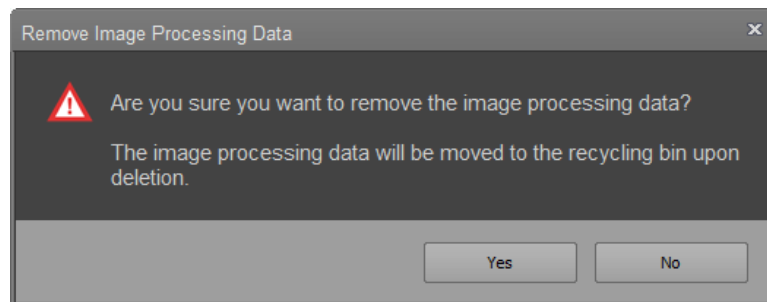
**IMPORTANT!** When image processing completes for each sample, any existing results are overwritten.

---

- If **Do this for all the samples** is selected and **Skip** is clicked, selected samples that already have image processing data are not added to the queue and only those samples that do not have existing data are processed.

## Remove Image Processing Data dialog

The **Remove Image Processing Data dialog** helps to prevent image processing data from being accidentally removed from Samples. The dialog is displayed when you select **Remove Image Processing Data** option from a **Sample context menu**.



- To remove the image processing data from the sample, click **Yes**. The image processing data are moved to the recycling bin.
- To close the dialog without removing the image processing data, click **No** or **X**.

## Link Account dialog

The **Link Account** dialog enables you to register the Attune™ Cytometric Software with the Connect cloud-based platform and link the instrument to the account on the **Instrument Connect** website.



---

**Note:** Only System Administrator and Administrator accounts can register the instrument with the Connect platform. After the device (instrument or software) is registered with the Connect platform, individual User accounts can link the registered instrument with their Connect accounts using the Cloud sign in feature. This allows a single instrument to be associated with multiple Connect accounts.

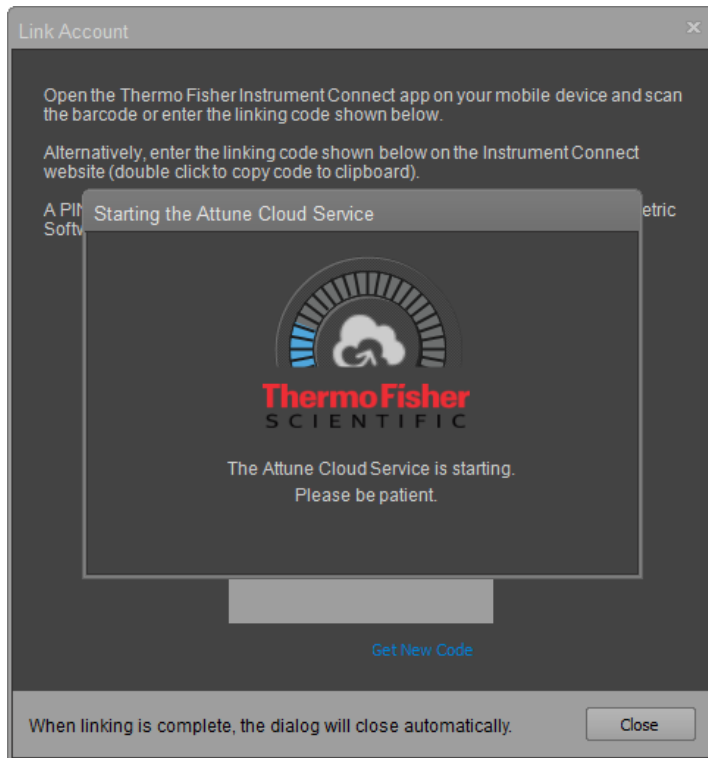
---

- The **Link Account** dialog provides two options to register the Attune™ Cytometric Software:
  - **Mobile:** Enables you to register the device software using a unique QR code with the QR code scanner in the **Instrument Connect** mobile application (“Link account with the QR code” on page 783).
  - **Linking Code:** Enables you to register the device software using a unique linking code from the Connect platform to register the device software on the **Connect website** (“Link account with the linking code” on page 785).
- If you click **Cancel** or the **X** button, the dialog closes without completing the device registration.

## Link account with the QR code

To register the device software with a unique QR code using the QR code scanner in the **Thermo Fisher Scientific Instrument Connect** mobile application:

1. On the **Options dialog** ▶ **Configuration tab** (“Configuration” on page 695), click **Register Device**. The software displays the **Starting the Attune™ Cloud Service™** dialog as the instrument establishes a connection.



2. When the connection with the cloud service is established, the **Link Account** dialog prompts you to link the cloud account by using either the **QR code** option (for mobile app devices) or a **linking code** (for web or mobile apps).

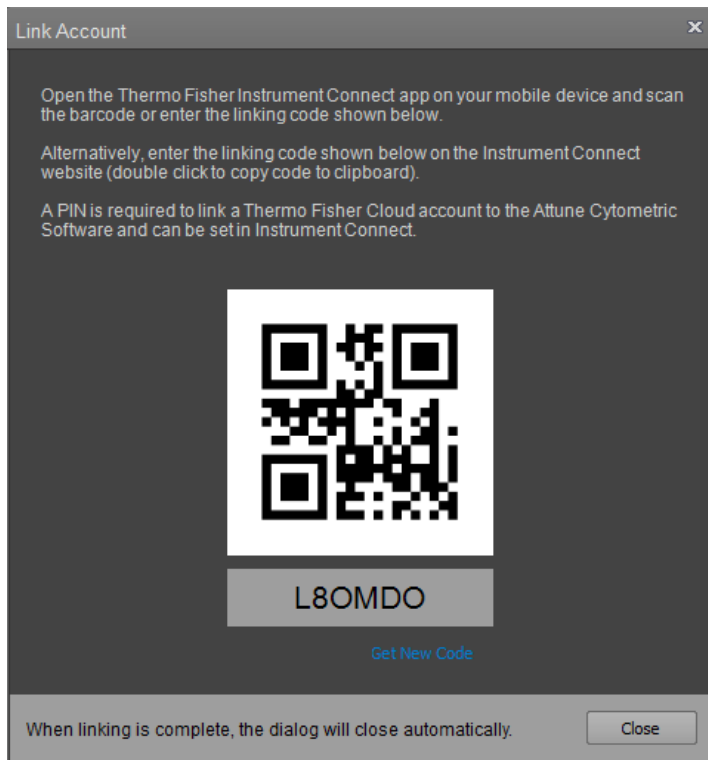


3. If an error occurs when retrieving the QR code, a warning banner is displayed on the dialog that provides a description of the error.  
To request another linking code, click the **Get New Code** hyperlink under the QR code. A new QR code is generated and displayed on the dialog.
4. When you have the QR code, open the **Thermo Fisher Scientific Instrument Connect** application on your mobile device, then scan the QR barcode image.
5. The **Instrument Connect** mobile application automatically sends the QR code to the Connect platform.
6. When registration is complete, the Attune™ Cytometric Software receives a successful response from the Connect platform, then the dialog automatically closes.

## Link account with the linking code

To register the device software using a unique linking code from the Connect platform:

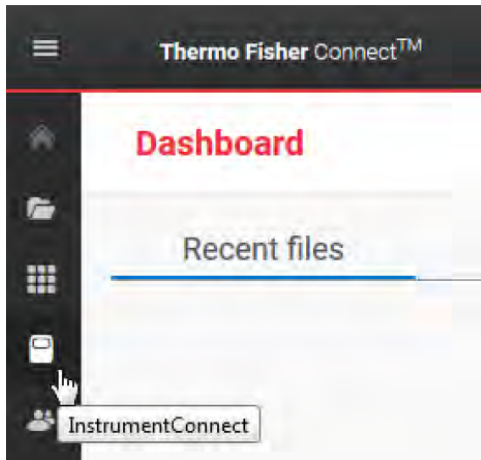
1. On the **Options dialog** ▶ **Configuration tab** (“Configuration” on page 695), click **Register Device**.
2. When the connection with the cloud service is established, double-click the **linking code** below the QR code to copy the linking code to the Windows™ clipboard.



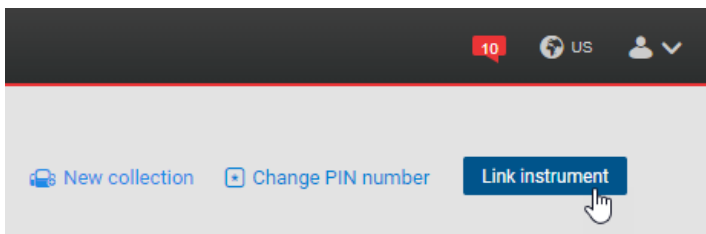
3. If an error occurs when retrieving the linking code, a warning banner is displayed on the dialog that provides a description of the error.

To request another linking code, click the **Get New Code** hyperlink under the linking code. A new linking code is generated and displayed on the dialog.

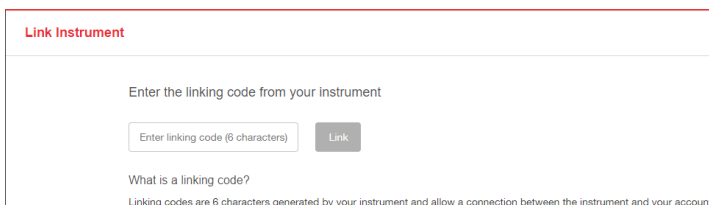
- When you have the linking code, go to [thermofisher.com](https://thermofisher.com), sign in to your Connect account, then click the **Instrument Connect** icon.



- In the Instrument tab, click **Link instrument** to open the **Link instrument dialog**.

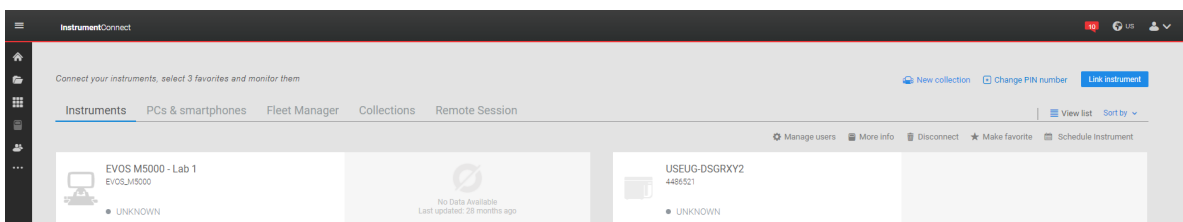


- Enter the unique linking code you have copied on the **Link Account dialog**, then click **Send**.



- When registration is complete, the Attune™ Cytometric Software receives a successful response from the Connect platform, then the dialog automatically closes.

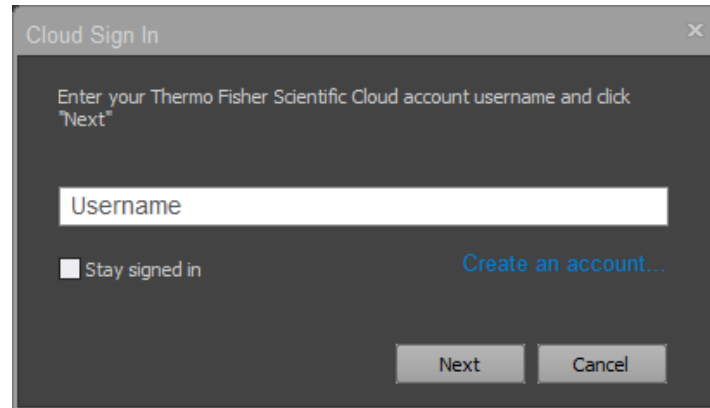
The registered device is listed in the Instruments section of your Connect account.





## Cloud Sign In dialog

The **Cloud Sign In** dialog enables you to sign in to your Connect account, Thermo Fisher Scientific's cloud-based platform, with your account username and password.

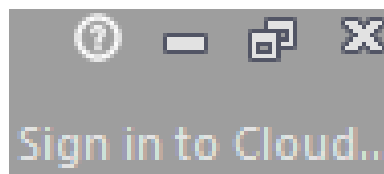


---

**Note:** Only System Administrator and Administrator accounts can register the instrument with the Connect platform. After the device (instrument or software) is registered with the Connect platform, separate User accounts can connect the registered instrument with their Connect accounts using the **Cloud Sign In**. This allows a single instrument to be associated with multiple Connect accounts.

---

1. To sign in to the Connect account, click **Sign in to Cloud** button, located at the top right corner of the screen. The **Cloud Sign In** dialog opens.

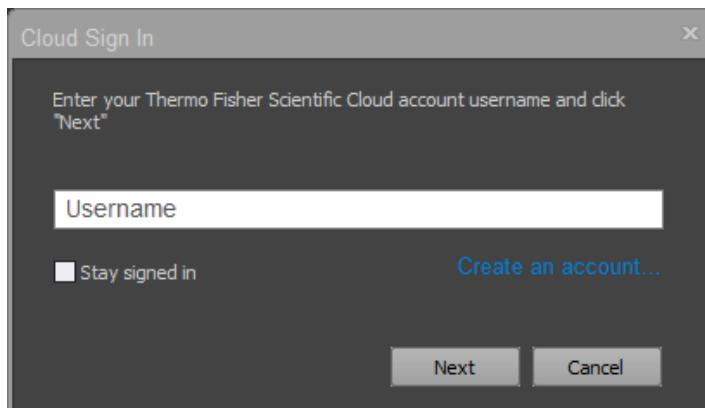


2. In the dialog, enter the Connect username in the **Username** field, then click **Next**.
  - If you do not yet have a Connect account, click **Create an account...** hyperlink, which opens a browser directed to the global registration site ([www.thermofisher.com/global-registration/registration](http://www.thermofisher.com/global-registration/registration)) to enable you to create an account.
  - To automatically sign in to the Connect platform the next time you sign in to the Attune™ Cytometric Software, select **Stay signed in**.  
If the **Stay signed in** checkbox is unchecked, every time you sign out of the Attune™ Cytometric Software, the sign in credentials are purged, which requires you to sign in the next time you want to connect to the Connect platform.

3. If you have not yet used your Connect account on the Attune™ Cytometric Software, the software displays the Link Account dialog (“Link Account dialog” on page 782), which enables you to connect the account to the software for the first time.

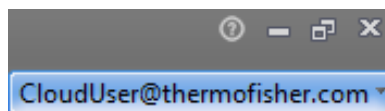


4. If you were the user who had registered the instrument or if you had already linked the instrument to the Connect account, the software prompts you to enter the **four-digit PIN** to connect the Connect account to the Attune™ Cytometric Software.

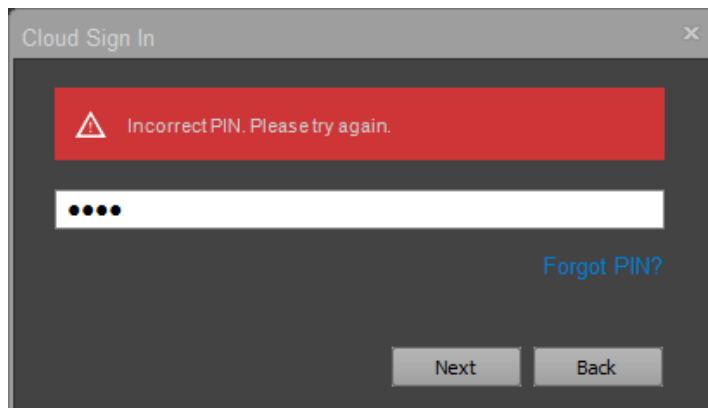


The **Forgot PIN?** hyperlink directs you to the web version of Instrument Connect application, where you can change or reset the PIN number.

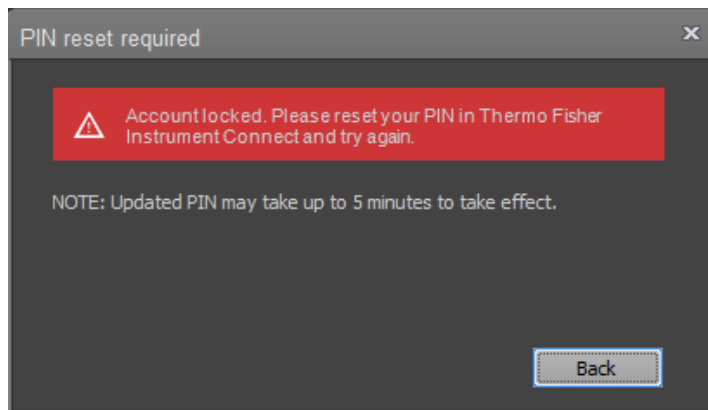
5. Providing a correct PIN number completes the account linking process and the **Sign in to Cloud** button located at the top right corner of the screen, becomes a dropdown button that displays the **Connect username** as the main button text.



6. If you provide a wrong PIN, you are notified and asked to try again.



7. If you provide the wrong PIN 5 times in a row, the account is locked and you will be required to reset the PIN in Instrument Connect before trying again.



After you enter a new PIN on the Thermo Fisher Scientific Instrument Connect website or mobile app, the **Incorrect PIN** dialog changes to the **Enter PIN** dialog.

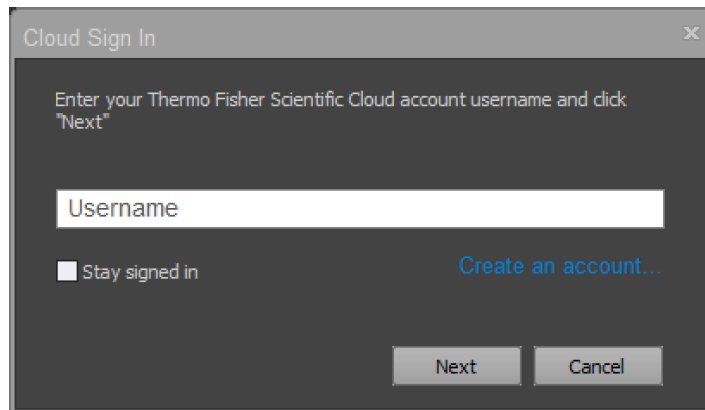
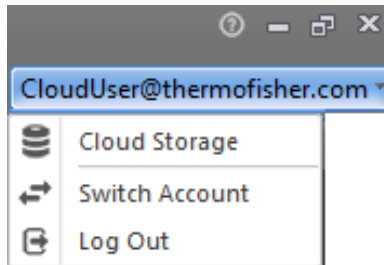
---

**Note:** The account lock out is only a local lock that prevents the user from linking their cloud account to the Attune™ Cytometric Software; it does not completely lock the Thermo Fisher Scientific Cloud account on the web.

---

## Switch Account dialog

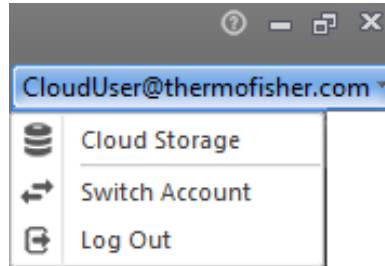
When you select **Switch Account** on the **Sign in to Cloud...** dropdown, the **Cloud Sign In** dialog (“Cloud Sign In dialog” on page 787) opens and enables you to sign in with a different username and for-digit PIN.



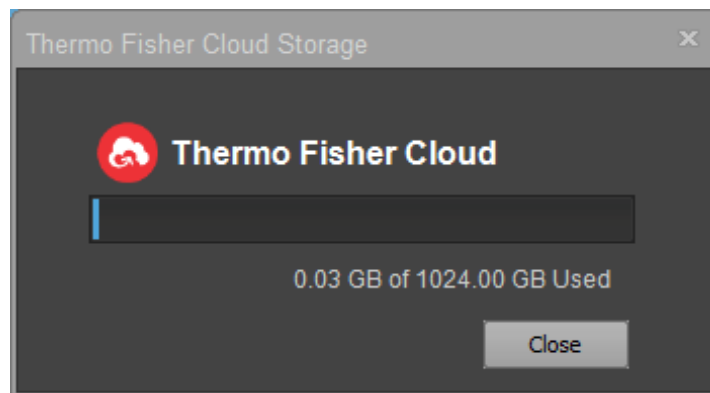
- The Switch account workflow is identical to that of Cloud Sign In, as described on “Cloud Sign In dialog” on page 787.
- When you switch to another Connect account successfully, the software automatically signs out of the old cloud account and clears the **Stay sign in** setting.

## Connect storage dialog

When you select **Cloud Storage** on the **Sign in to Cloud...** dropdown list, the **Connect Storage** dialog opens and shows the available storage on the Connect account for the current user.

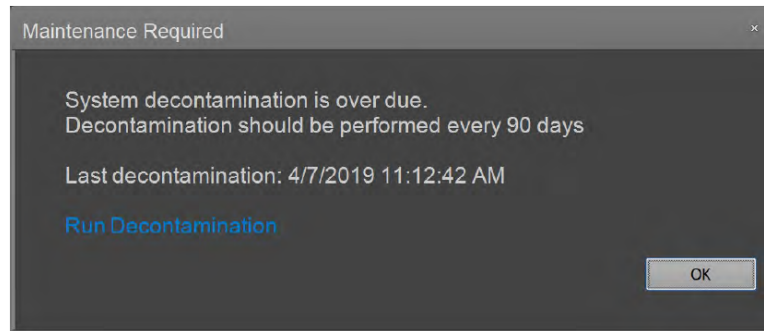


The progress bar on the dialog graphically displays the total used space.



## Maintenance Required dialog

If there is overdue maintenance and the instrument is online and connected, the **Maintenance Required** dialog is displayed during the sign in and describes which maintenance procedure is overdue.

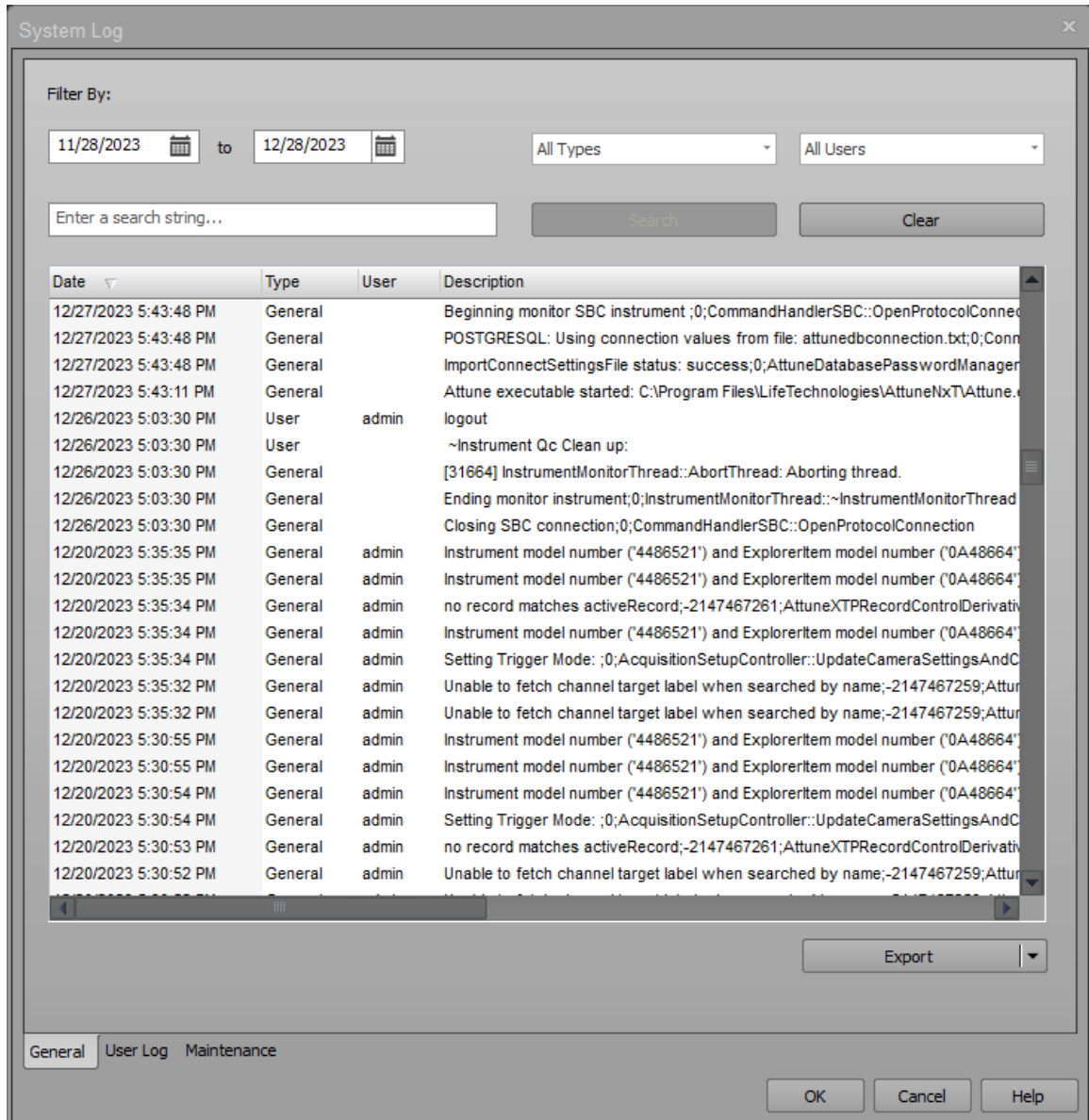


- To start the described maintenance procedure and close the dialog, click the **maintenance procedure hyperlink** that is displayed in the dialog. The hyperlink is displayed only for users who have permission to perform the maintenance procedure.
- To dismiss the dialog without performing the maintenance procedure, click **OK**. However, the dialog will be displayed at each subsequent sign in until the required maintenance procedure is performed.

## System Log dialog

**System Log dialog** is displayed when the **System Log** or the **Maintenance Log** button is clicked on the **Instrument** ribbon tab (see "Instrument tab" > "Functions group" on page 90).

- Click **System Log** in the **Instrument** ribbon tab to open the dialog to the **General** tab.
- Click **Maintenance Log** in the **Instrument** ribbon tab to open the dialog to the **Maintenance** tab.



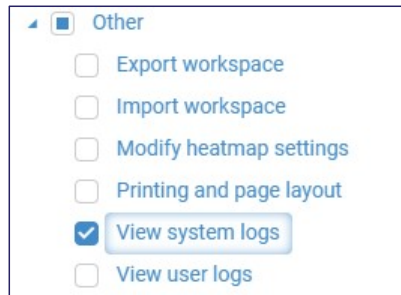
The **System Log dialog** enables users to view the system transactions and maintenance history based on the user account privileges described in "Account types" ("Default local Attune™ accounts" on page 42).

Depending on the account type, the **System Log dialog** displays different information:

- For **Service** accounts, the **User column** and **User combo box** in the **General** tab, and the **User Log** tab are not visible.
- For **System Administrator** and **Administrator** accounts, the **Service column** in the **General** tab is not visible. Both **System Administrator** and **Administrator** accounts have access to the **User Log** and can view all user level information.
- For general **user** accounts, the **Service column**, **User column**, and **User filter** in the **General** tab, and the **User Log** tab are not visible.
- All user accounts can see and access the **Maintenance** tab.
- Only **System Administrator**, local **Administrator**, and **SAE Users** with the permission to **Configure Security and Auditing** can view the entirety of the **User column** and sort by **All** users. Otherwise, they can only select themselves as an option and see other users as a representative GUID for privacy.

---

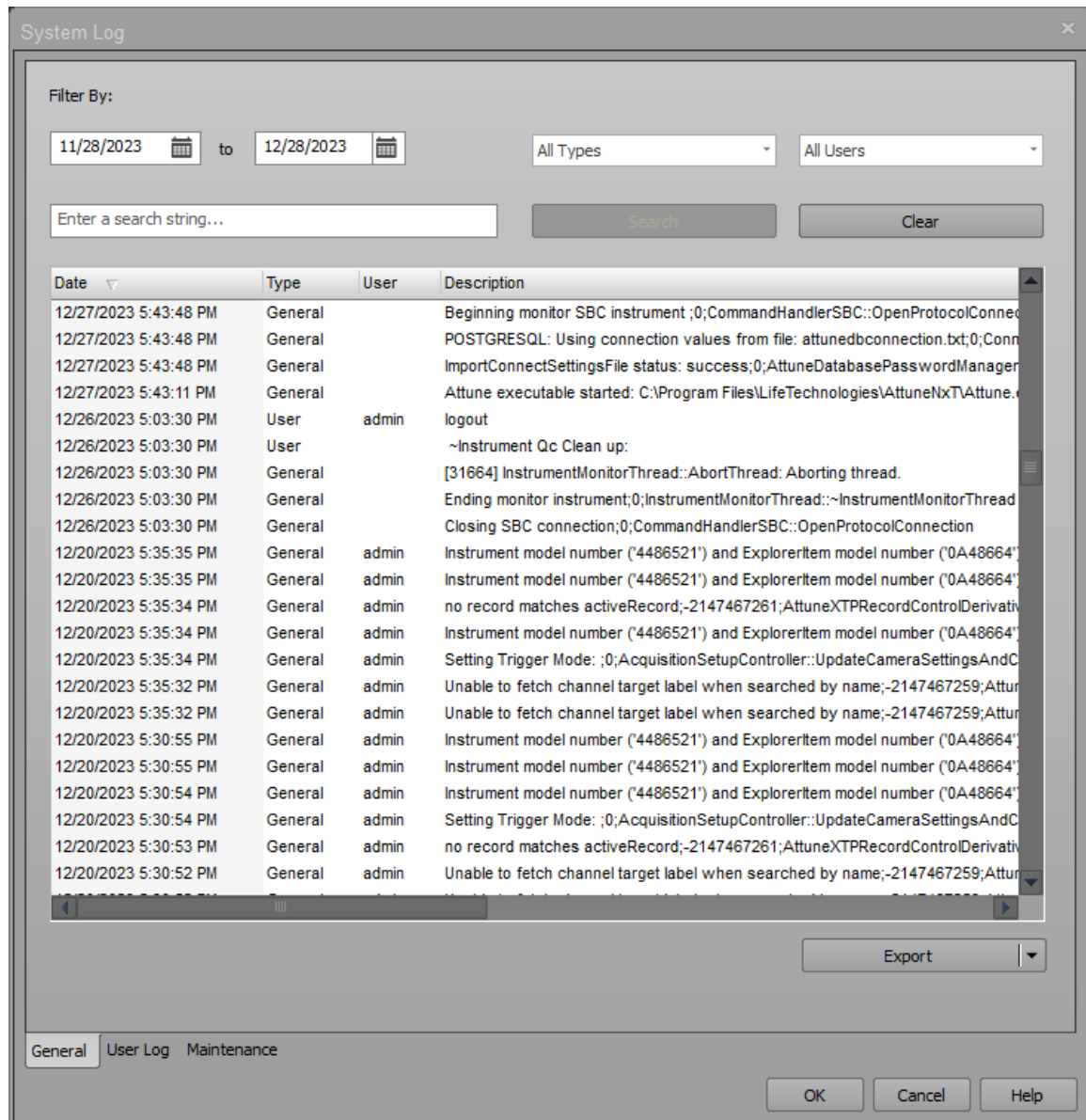
**Note:** In the **SAE mode**, you need to have **Attune Software > Other > View system logs** permission selected to view the **Maintenance** tab (see “SAE account permissions” on page 887).





## General tab

**General** tab of the **System Log** dialog enables you to view general system transactions based on the account privileges described in “Account types” (“Default local Attune™ accounts” on page 42). It contains the **filters**, the **log table**, and the **Export/Print** split button.



**Note:** Only **System Administrator**, local **Administrator**, and **SAE Users** with the permission to **Configure Security and Auditing** can view the entirety of the **User** column and sort by **All** users. Otherwise, they can only select themselves as an option and see other users as a representative GUID for privacy.

## Filters

Filters are available based on account type. For all filter types, the filter is applied to the list after the filter choice is made. The filters reset when **Clear** is clicked.

- **Date range:** Selects the date range for the system log entries.
- **Type:** Contains the **Error**, **Service**, **Info**, **User**, and **All** options. By default, **All** is selected.
- **User:** Contains all users of the system. Users are as their full name (last, first), and the options include **All Users**. By default, **All Users** is selected.
- **Search box** and **Search button:** **Search box** enables you to search through any column to find rows containing the phrase typed in the box. Clicking the **Search** button applies the search prompt to the list. If no text is entered in the search textbox, the **Search** button is disabled.
- **Clear:** Clears all filters and resets the filters to their defaults including the entered search text.

## Log table

The columns within the Log table depend on the user type. At a minimum, the following columns exist: **Date**, **User**, **Type**, **Description**, and **Service**.

The log table contains the following information, depending on the user type.

Action	Type	Description	Service Value
Performance Test	Info	Result	N/A
Baseline	Info	Result	N/A
Functions (wash, rinse, decontaminate, etc.)	Info	Function and function cycle	N/A
Plate run time	Info	<Plate experiment name> HH:MM:SS	N/A
Sample run time	Info	<sample name> HH:MM:SS	N/A
Error messages that need customer intervention (low fluid level, etc.)	Info	Which Fluid low warning triggered	N/A
Error messages that generate an error dialog and error number	Error	Error number	The error description as retrieved from instrument API
Log in	User	Username login	N/A

(continued)

Action	Type	Description	Service Value
Log out	User	Username logout	N/A
PZT calibration	Service	PZT calibration performed and outcome	Message: Resonance Power Value or Error message
PZT Acoustic Power Adjustment	Service	PZT acoustic power adjusted	Flow Rate: Value
Laser Delay Adjustment	Service	Laser Delay Adjusted for <color> laser	Value
System Test	Service	Outcome	Specific errors

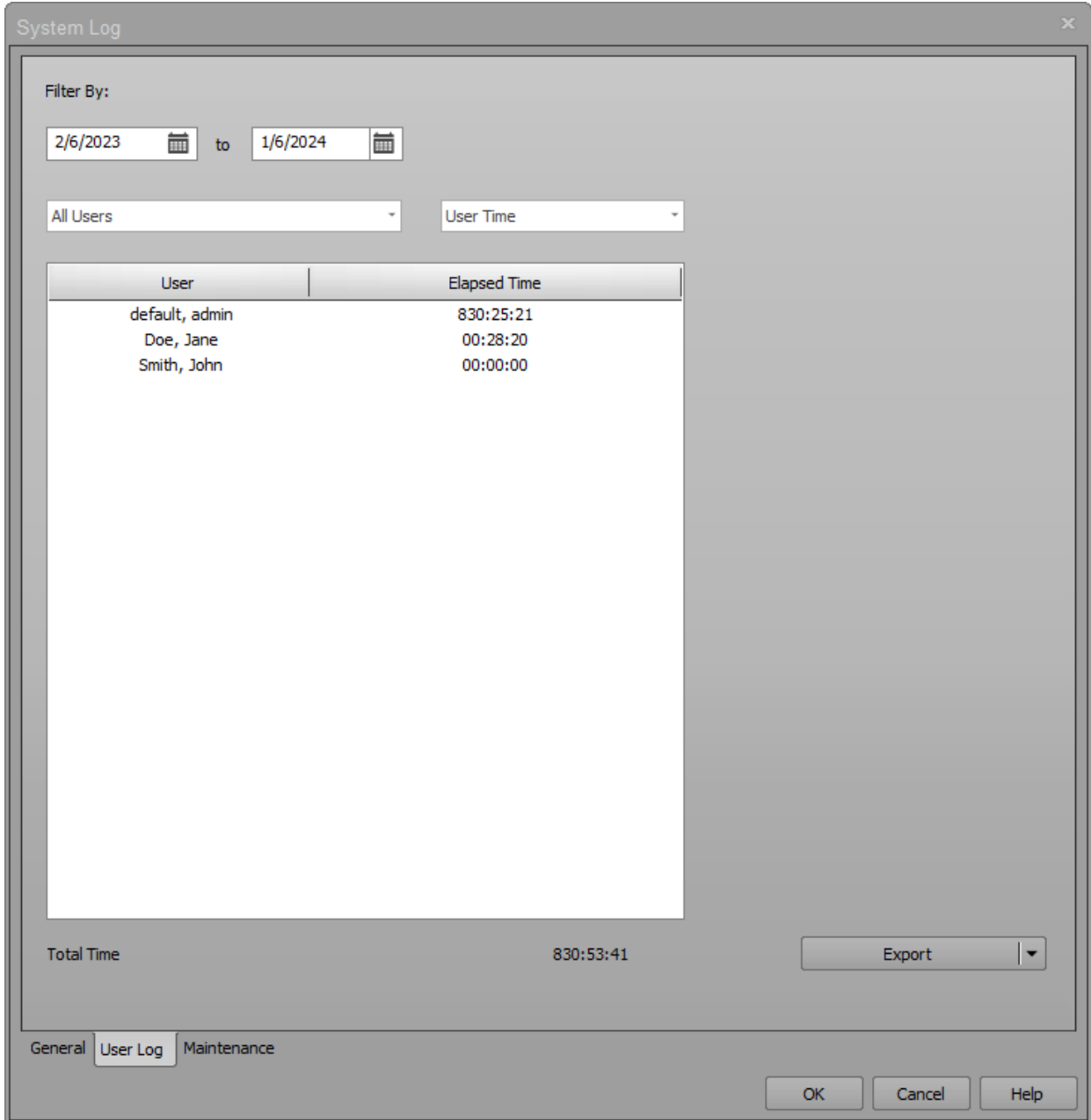
### Export/Print split button

The **Export/Print** split button enables you to export or print the **System log**.

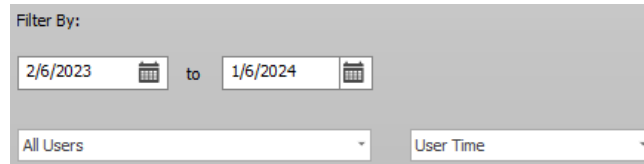
- Clicking the **Export** part of the button opens the **File Save (Export)** dialog (“File Save (Export) dialog” on page 715). The exported file is saved as an ALF file.
- The dropdown arrow of the split button enables you to select **Print**, which opens the **Print** dialog (“Print dialog” on page 730).

## User Log tab

**User Log** tab of the **System Log dialog** shows the elapsed login time for any user within a specified date range.



## Filters



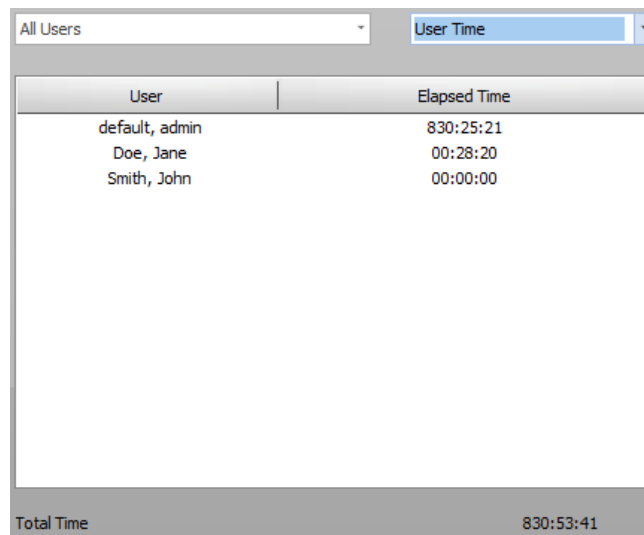
Filter By:

2/6/2023 to 1/6/2024

All Users User Time

- **Date range:** Selects the date range for the log entries.
- **User:** Contains the list of all users of the system in a dropdown. Users are listed with their full name (last, first), and the dropdown list also includes **All Users**. By default, **All Users** is selected.
- **User Time:** Contains the options to show elapsed **User Time** or **Sample Count** for each user listed in the **Log Table**. The default option is set to **User Time**.

## Log table



User	Elapsed Time
default, admin	830:25:21
Doe, Jane	00:28:20
Smith, John	00:00:00
Total Time	830:53:41

The table displays columns according to the choice in the **User Time** dropdown:

- When **User Time** is selected, the columns are **User** and **Elapsed Time**.
- When **Sample Count** is selected, the columns are **User** and **Sample counts**.

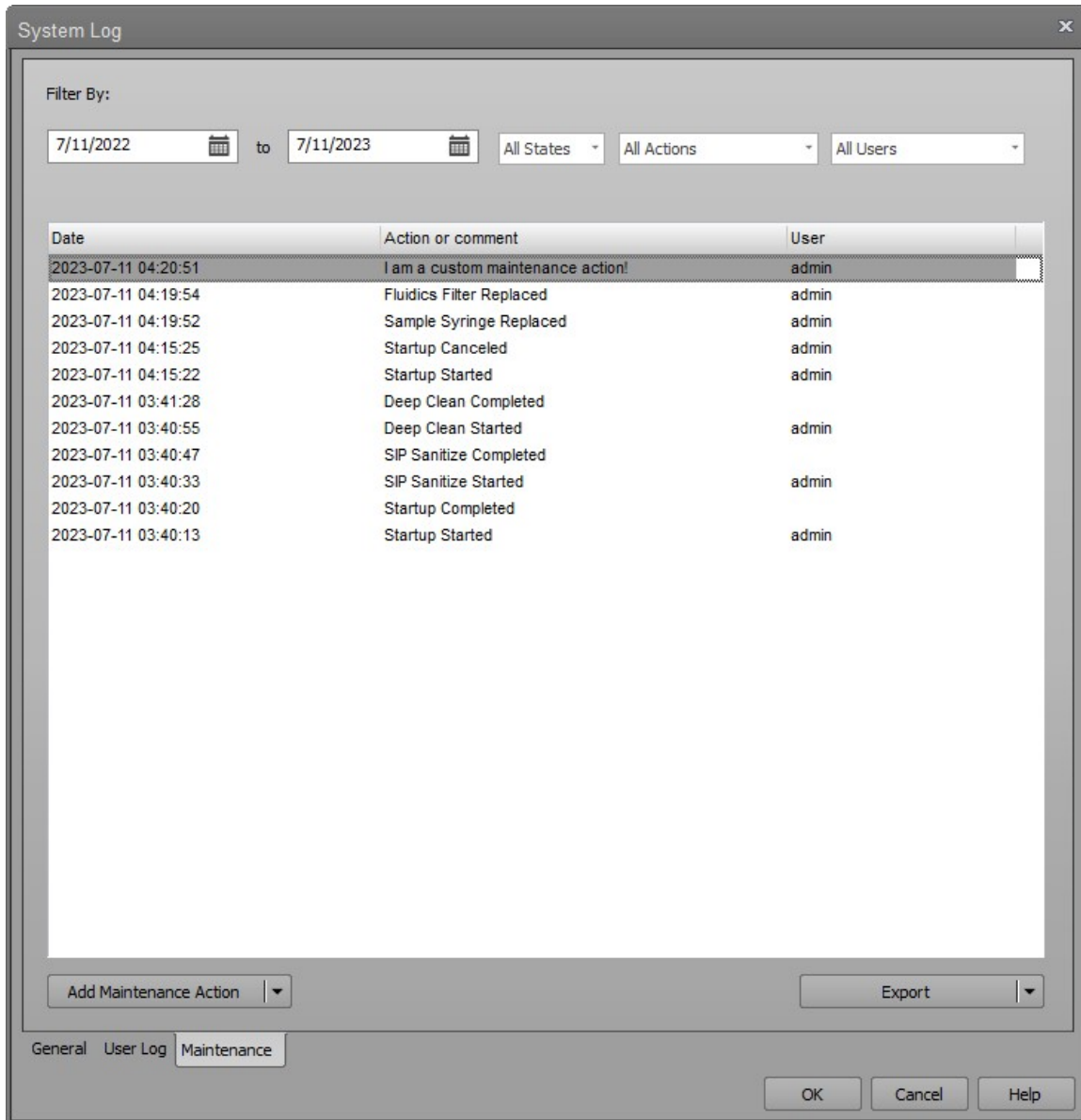
## Export/Print split button

The **Export/Print** split button enables you to export or print the **User log**.

- Clicking the **Export** part of the button opens the **File Save (Export)** dialog (“File Save (Export) dialog” on page 715).
- The dropdown arrow of the split button enables you to select **Print**, which opens the **Print** dialog (“Print dialog” on page 730).

## Maintenance tab

**Maintenance** tab of the **System Log dialog** provides a record of the instrument maintenance tasks performed by all users and when they were performed. It also enables you to add custom maintenance tasks that are not automatically included to the **Maintenance Viewer**. The **Maintenance Viewer** can be sorted by **Date**, **Action or comment**, and **User**, and its contents can be exported locally or printed to keep record of the instrument maintenance history.

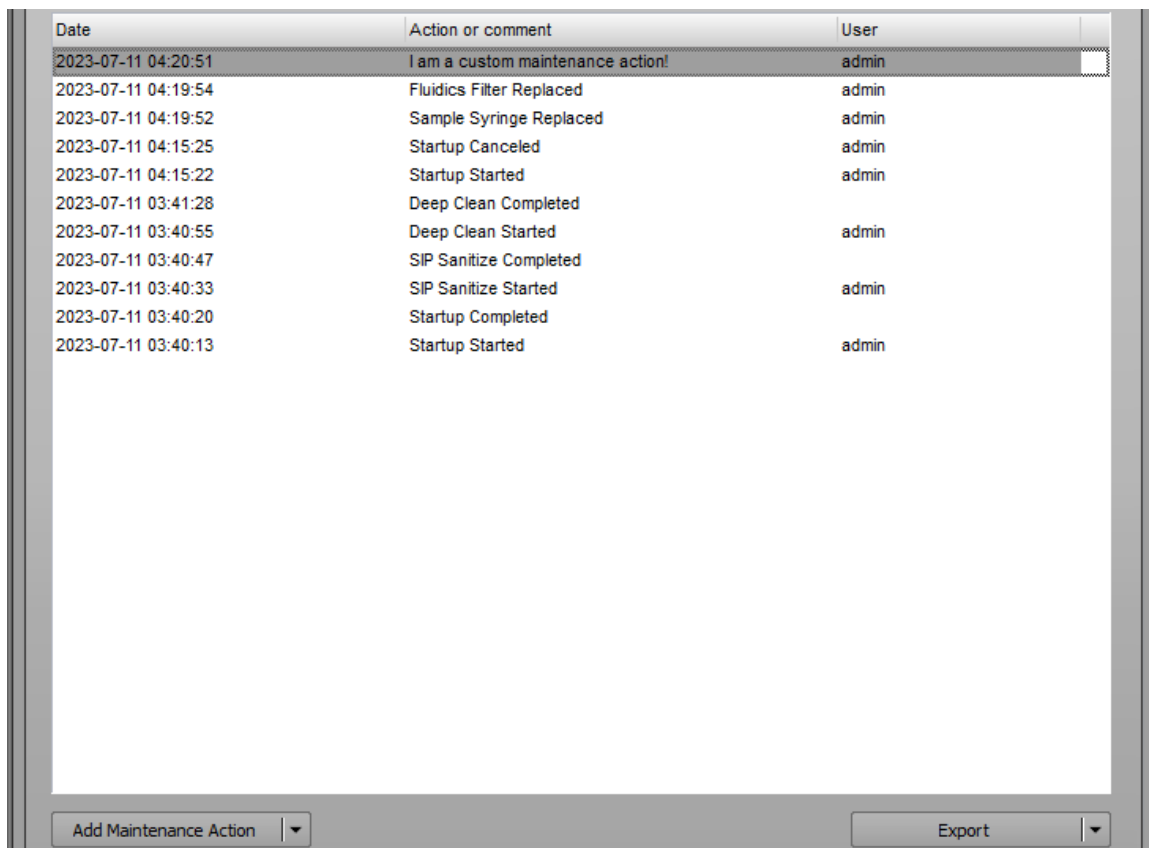


**Note:** Only local **Administrator**, **System Administrator**, and **SAE Users** with the permission to **Configure Security and Auditing** can view the **User** column and sort by **User**.

**Note:** In the **SAE mode**, you need to have **Attune Software** ▶ **Other** ▶ **View system logs** permission selected to view the **Maintenance** tab and to add custom maintenance tasks (see “SAE account permissions” on page 887).

## Maintenance Viewer

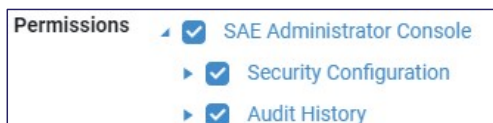
The **Maintenance Viewer** lists the instrument maintenance tasks that were performed, which can be sorted by **Date**, **Action or comment**, and **User**. The list can also be exported locally or printed to keep record of the instrument maintenance history.



Date	Action or comment	User
2023-07-11 04:20:51	I am a custom maintenance action!	admin
2023-07-11 04:19:54	Fluidics Filter Replaced	admin
2023-07-11 04:19:52	Sample Syringe Replaced	admin
2023-07-11 04:15:25	Startup Canceled	admin
2023-07-11 04:15:22	Startup Started	admin
2023-07-11 03:41:28	Deep Clean Completed	
2023-07-11 03:40:55	Deep Clean Started	admin
2023-07-11 03:40:47	SIP Sanitize Completed	
2023-07-11 03:40:33	SIP Sanitize Started	admin
2023-07-11 03:40:20	Startup Completed	
2023-07-11 03:40:13	Startup Started	admin

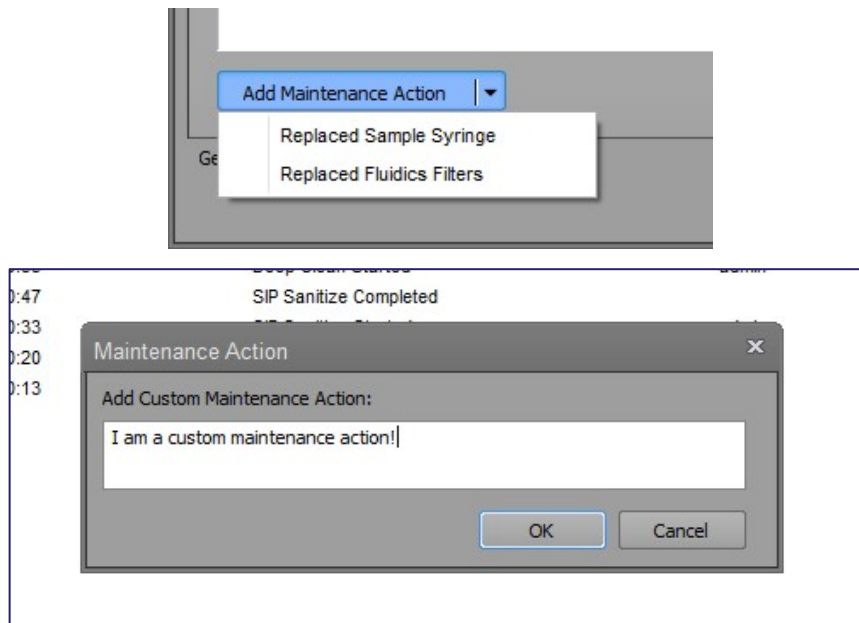
At the bottom of the dialog, there is an "Add Maintenance Action" button on the left and an "Export" button on the right.

**Note:** Both the **Security Configuration** and the **Audit History** permissions must be selected in the **SAE Administrator Console** (see “SAE account permissions” on page 887) to see the **User** column in the **Maintenance Viewer**, and to view which user has performed which instrument function.



- The **Maintenance Viewer** automatically logs the following maintenance tasks when they are performed:
  - Startup
  - Shutdown
  - Decontamination
  - SIP Sanitize (Tube and Autosampler)
  - Deep Clean
  - Debubble
  - Unclog
  - Autosampler Calibration
  - PZT Calibration

- The **Maintenance Log** does not record the following maintenance tasks, but they can be manually entered into the log (see Add Maintenance Action).
  - Rinse
  - Test Plate
  - Self Test
  - Run/Record of Sample
  - Sample Recovery
  - Service Camera
  - Gear Pump Calibrations
- When an instrument maintenance function is started, the **Maintenance Viewer** records the timestamp when it starts, the user, and the action.
- When the function is complete, the **Maintenance Viewer** records the timestamp and notes the completion of the same function. The user is not recorded to minimize redundancy in the log.
- When an instrument function is canceled by selecting **Stop** in the **Instrument** ribbon tab, the **Maintenance Viewer** records the timestamp of cancellation, the user, and the function that was canceled.
- The **Maintenance Viewer** does not record instrument functions that are stopped when an error occurs.
- **Add Maintenance Action** enables you to log maintenance actions that are not recorded automatically. You can select the action from the dropdown list or enter a custom maintenance action directly.



---

**Note:** When adding a custom maintenance action, there is a 100-character limit. The characters used for the custom maintenance action cannot include an ampersand (&).

---



## Filters

Filters enable you to filter instrument maintenance record by **Date range**, **State**, **Action**, and **User**. For all filter types, the filter is applied to the list after the filter choice is made.



The screenshot shows a 'Filter By:' section with three main components: a date range selector, a state dropdown, an action dropdown, and a user dropdown. The date range is set from 7/11/2022 to 7/11/2023. The state dropdown is set to 'All States', the action dropdown is set to 'All Actions', and the user dropdown is set to 'All Users'.

- **Date range:** Selects the date range for the system log entries.
- **State:** Contains the **Error**, **Service**, **Info**, **User**, and **All States** options. By default, **All States** is selected.
- **Action:** Contains all actions that are listed in the **Maintenance Viewer**. By default, **All Actions** is selected.
- **Users:** Contains all users of the system, including SAE users. Users are listed by their username and the dropdown choices also include **All Users**. By default, **All Users** is selected. In the SAE mode, the users are listed by their SAE usernames, if correct permissions allow their viewing.

## Export/Print split button

The **Export/Print** split button enables you to export or print the **Maintenance log**.

- Clicking the **Export** part of the button opens the **File Save (Export)** dialog (“File Save (Export) dialog” on page 715).
- The dropdown arrow of the split button enables you to select **Print**, which opens the **Print** dialog (“Print dialog” on page 730).

# SAE Signing dialogs

## Overview

**SAE Signing dialogs** include the dialogs listed below, which are displayed when the corresponding **SAE Signing** button is clicked on the **SAE** ribbon tab (“SAE tab” on page 114):

**Request Signature** (“Request Signature dialog” on page 806)

**e-Signature Request** (“e-Signature Request dialog” on page 807)

**Pending Signatures** (“Sign Records dialog” on page 808)

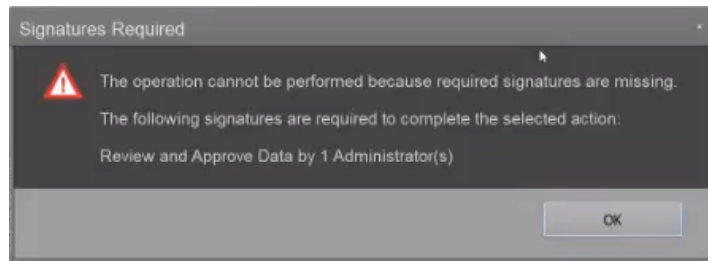
**Sign Record** (“Sign Experiment dialog” on page 810)

**e-Signature Record Report** (“e-Signature Record Report History dialog” on page 812) dialogs, which are .

You can configure the **e-Signature settings** (actions requiring signatures, number of signatures required for each action, reasons available for e-Signature, data to be signed) in the **e-Signature** tab of the **SAE Administrator Console** (“e-Signature tab” on page 909).

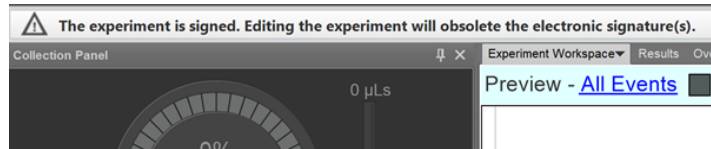
## e-Signature workflow

- Actions that require signatures to be completed cannot be performed unless the required signatures are provided.
- **Request Signature** dialog (“Request Signature dialog” on page 806) displays the **Report** of the active Experiment and enables you to request signatures for the selected action. Until it is signed, the **Report** displays the watermark that states **Unsigned**.
- If there are pending e-Signature requests, the **e-Signature Request** dialog (“e-Signature Request dialog” on page 807) is displayed on signing into the software. The **e-Signature Request** dialog also contains a link to the **Sign Records** dialog.
- If you attempt to perform an action that has unmet signature requirements, the **e-Signatures Required** dialog is displayed. The dialog shows the number outstanding roles and meanings that must be signed for before you can perform the action.



- The **Sign Records** dialog (“Sign Records dialog” on page 808) enables you to view and sign pending signature requests. After the report has been signed, the **Experiment Report** removes the watermark and changes the report status to **Current**.

When all signatures are current, the notification bar on the Attune™ desktop states **The experiment is signed. Editing the experiment will obsolete the electronic signature(s).**

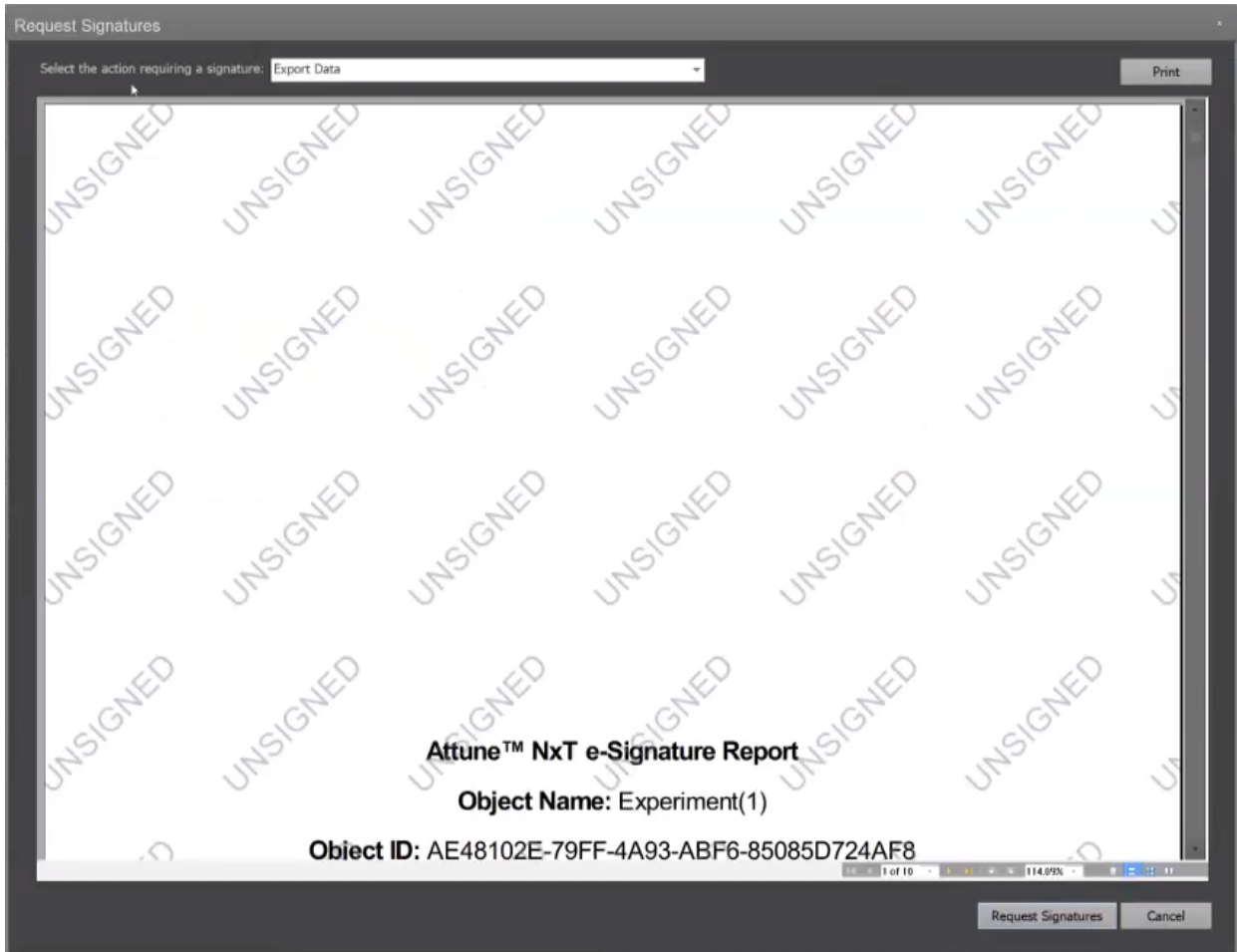


- Signatures represent the state of the Experiment at the time of signing. If any modifications are made to the Experiment after signing, the signatures become obsolete and the **Experiment Report** displays the watermark that states **Obsolete**. In such cases, actions that require signatures cannot be performed unless new signatures are provided.
- **e-Signature Record Report History** dialog (“e-Signature Record Report History dialog” on page 812) lets you view and print signed records.

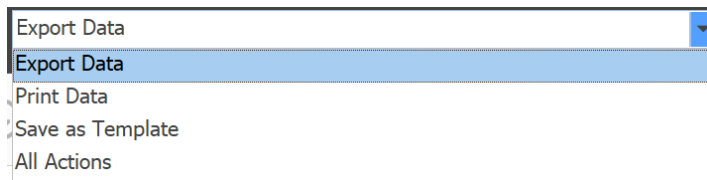
## Request Signature dialog

**Request Signature** dialog displays the **Report** of the active Experiment and enables you to request signatures for the current experiment.

To view this dialog, click the **Request Signatures** button on the **SAE** Ribbon tab (“SAE tab” on page 114).



1. Select the desired action from the **Select the action requiring a signature** dropdown menu.



---

**Note:** Only actions that have been configured to require signatures are displayed in the **Select the action requiring a signature** dropdown menu (see “Select the actions that require e-Signature” on page 910).

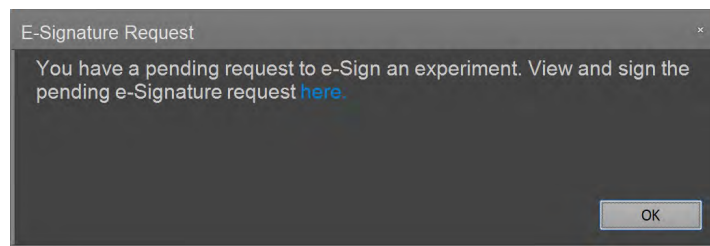
---

2. (Optional) Click **Print** to print the report for the Experiment.
3. Click **Request Signatures** to send an e-Signature request.

## e-Signature Request dialog

**e-Signature Request** dialog is displayed on signing into the software if there is a pending signature request needed of the user signing in.

The dialog is displayed each time a user signs in as long as there is a pending signature request.

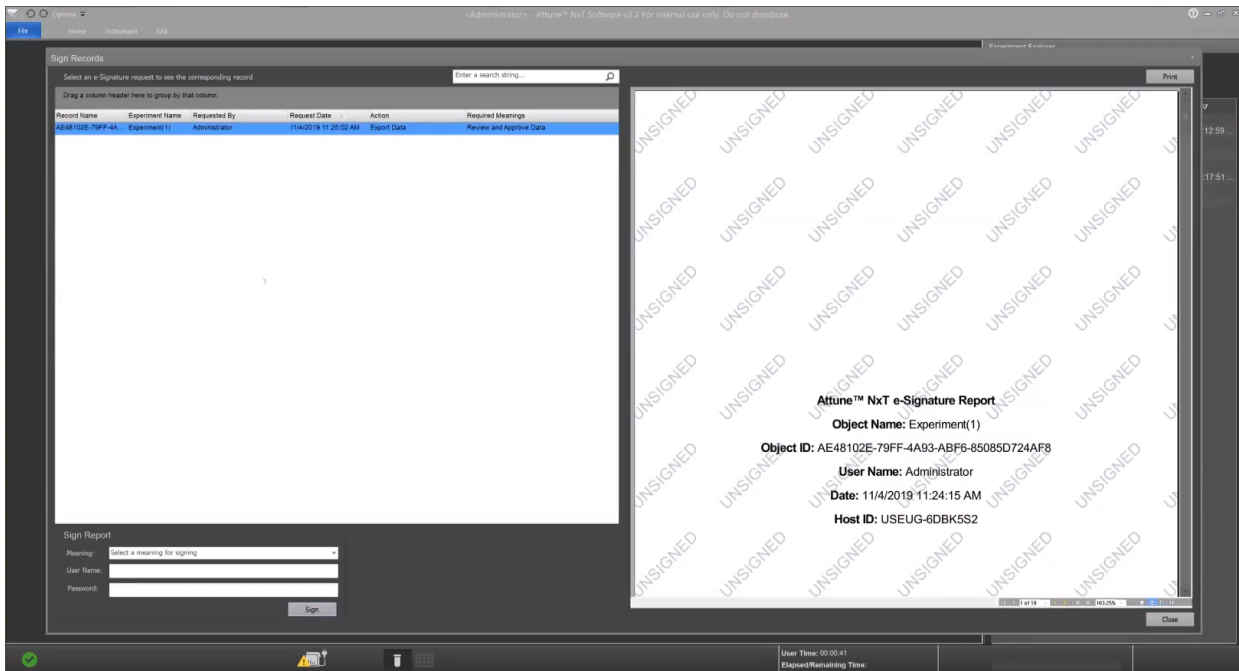
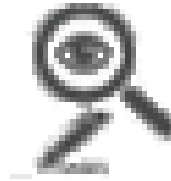


- Click the **here** hyperlink to close the dialog and open the **Sign Records** dialog (“Sign Records dialog” on page 808).
- Click **OK** to close the dialog without taking any further action.

## Sign Records dialog

**Sign Records** dialog enables you to view and sign pending signature requests.

To view the **Sign Records** dialog, click the **View Pending Signatures** button on the **SAE Ribbon** tab (“SAE tab” on page 114).



- The pending signature request is shown in the **Pending Signatures** panel on the left, which lists the Record name, Experiment name, the requester, the request date, requested action, and required meanings for the requested signature.

Record Name	Experiment Name	Requested By	Request Date	Action	Required Meanings
CBE14949-76BA-4...	Experiment(1)	admin	3/20/2019 5:28:32 PM	Export Data	Review and Approved Data   Review and Approv...

1. Select the record you want to sign for from the **Pending Signatures** panel.  
The **Experiment Report** for the selected signature request is displayed on the right panel.
2. (Optional) Click **Print** to print the report for the experiment.

3. Select the **Meaning** for the signature request in the **Sign Report** panel (which is listed in the **Required Meanings** column of the **Pending Signatures** panel), then enter the **Username** and **Password**.



The image shows a dialog box titled "Sign Report". It contains three input fields: a dropdown menu labeled "Meaning:" with the text "Select a meaning for signing" and a downward arrow; a text input field labeled "User Name:"; and another text input field labeled "Password:".

4. Click **Sign** to provide the e-Signature for the selected signature request.

## Sign Experiment dialog

**Sign Experiment** dialog shows the **Report** of the Experiment for which the signature is requested and enables you to sign the Experiment.

To view the **Sign Experiment** dialog, click the **Sign Record** button on the **SAE Ribbon** tab (“SAE tab” on page 114).



Sign Experiment(1)

Attune™ NxT Software Page 3

**Workspaces -- Group | 200 ul min | Experiment-level WS --- Page 1 of 1**

All Events - 200 ul min All Events - 200 ul min

Experiment: Experiment(1)  
Group: Group  
Sample: 200 ul min  
Time Recorded: 00:31:22

Sign Report

Meaning:

User Name:

Password:



1. To sign the **Report** for the Experiment, select the **Meaning** for the signature, then enter the **Username** and **Password**.



Sign Report

Meaning:

User Name:

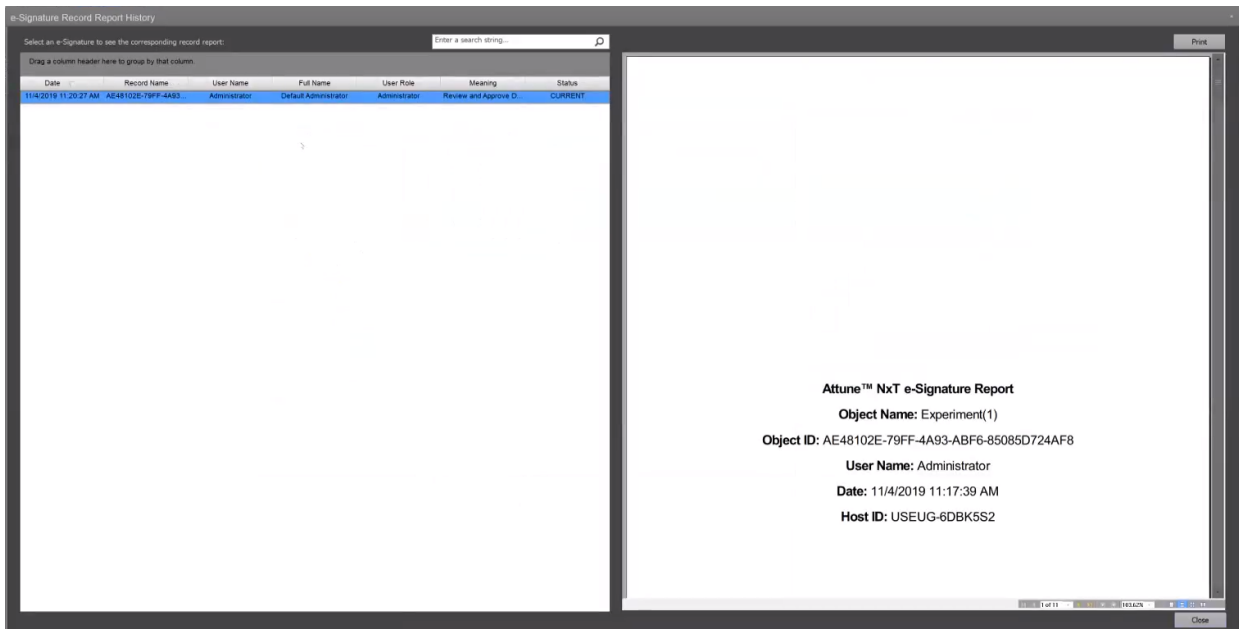
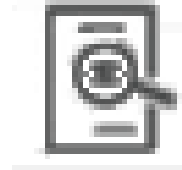
Password:

2. (Optional) Click **Print** to print the **Report** for the Experiment.
3. Click **Sign** to provide the e-Signature for the active Experiment Report.

## e-Signature Record Report History dialog

**e-Signature Record Report History** dialog enables you to view and print signed records.

To view the **e-Signature Record Report History** dialog, click the **View Signed Records** button on the **SAE Ribbon** tab (“SAE tab” on page 114).



- The e-Signatures that are on record are shown in the **Signatures** panel on the left, which contains the **Date**, **Record name**, **Username**, **Full name**, **User role**, **Meaning**, and **Status** columns.
- To search for an e-Signature record, enter a search string into the **search** field, then click **Enter**.
- To sort e-Signature records, drag the **column header** to the area where it states **Drag a column header here to group by that column**.
- To print an e-Signature report, select the report from the **Signatures** panel on the left, then click **Print** to print the report.

## Overview

The Attune™ Database Utility is a data backup program that is provided with the Attune™ Cytometric Software. This utility program enables you to back up the data to a custom location, either on-demand or at a specified, scheduled frequency.

The Attune™ Database Utility is used for the following procedures:

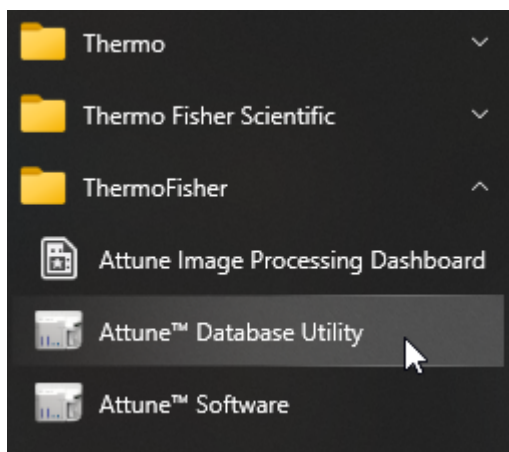
- **Backup** existing data and the database.
- **Restore** backed up data and database.
- **Schedule** an automate backup of the data and database.
- **Reinstall** the database.

## Startup

### Main application startup

The Attune™ Database Utility can be launched via the Windows™ **Start** menu or directly from the Attune™ Cytometric Software.

1. Select **Start** ▶ **All Programs** ▶ **ThermoFisher** ▶ **Attune™ Database Utility**.



Alternatively, start the Attune™ Database Utility from in the Attune™ Cytometric Software using the **Options** ▶ **Administrator options** ▶ **Backup options** dialog as described on “Data Management Options” on page 677.

2. The application opens to the **Login screen** (“Login screen” on page 814).

## Startup errors

### PostgreSQL error

If the database application is missing (PostgreSQL), the utility will not start and will display an error dialog with the following message:

“PostgreSQL 9.3.5 was not found. Please install the PostgreSQL using the Attune™ installer.”

### Database missing

If the Attune™ database is not present when starting the Attune™ Database Utility, the application launches in Recovery™ mode.

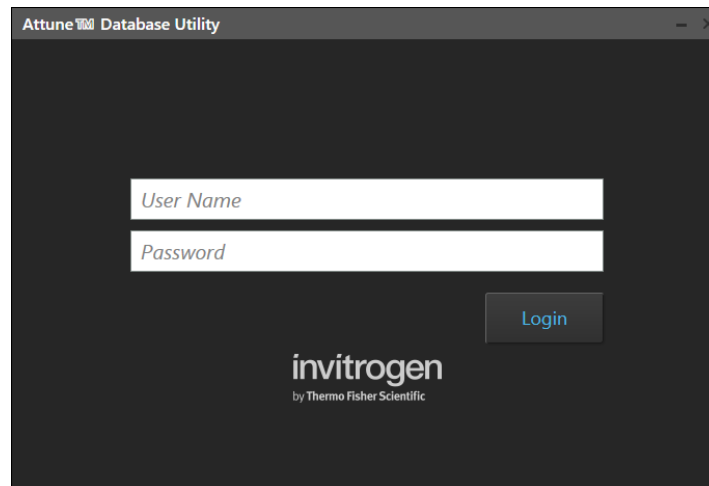
In Recovery mode, the application launches to the Restore™ page, which allows you to run the Restore™ or Reinstall database options. In this case, no login is required.

## Login screen

### Overview

The **Attune™ Database Utility Login** screen is the first screen that is displayed after the start of the utility. Each user must sign in before being allowed to use the utility.

Depending on the type of user (i.e., **Administrator**, **System Administrator**, **User**, etc.), specific features are available in the utility.




## User sign in

1. To sign into the **Attune™ Database Utility**, type a valid username and password in the corresponding field, then click **Login**.

Alternatively, press the **Enter** key on the computer keyboard after typing the username and password.

2. If the correct username and password are entered, the user login is completed and the **Attune™ Database Utility** opens.


If an invalid username or password is entered, the warning banner displays **Invalid username or password** (see “Login Warnings”).



△ Invalid username or password.

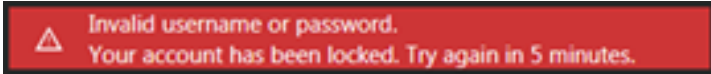
## Login warnings

- If an invalid username or password is entered, the warning banner displays **Invalid username or password**. A valid username and password must be entered to proceed.



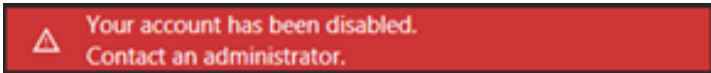
△ Invalid username or password.

- If the security settings (“System Security tab” on page 692) have been set to lockout a user after a specified number of failed attempts, the warning banner displays **Invalid username or password. Your account has been locked. Try again in X minutes.**



△ Invalid username or password.  
Your account has been locked. Try again in 5 minutes.

- If an attempt is made to login to an account that has been disabled, the warning banner displays **Your account has been disabled. Contact an administrator.**



△ Your account has been disabled.  
Contact an administrator.

# Main application

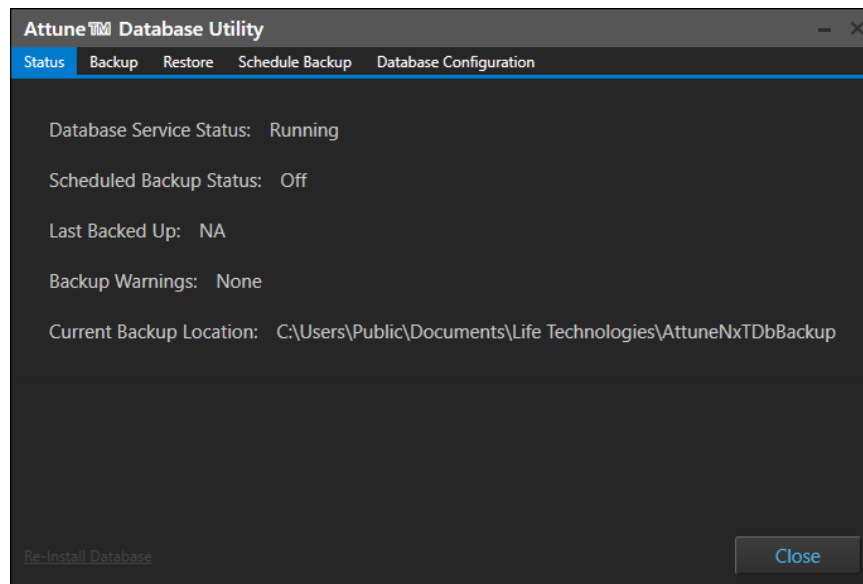
## Overview

The main application of the Attune™ Database Utility consists of a dialog window that contains the following controls:

- **Status** tab (“Status tab” on page 817)
- **Backup** tab (“Backup tab” on page 818)
- **Restore** tab (“Restore tab” on page 820)
- **Schedule Backup** tab (“Schedule Backup tab” on page 822)
- **Reinstall Database** hyperlink (“Reinstall Database” on page 823)
- **Database Configuration** tab (“Database Configuration tab” on page 824)

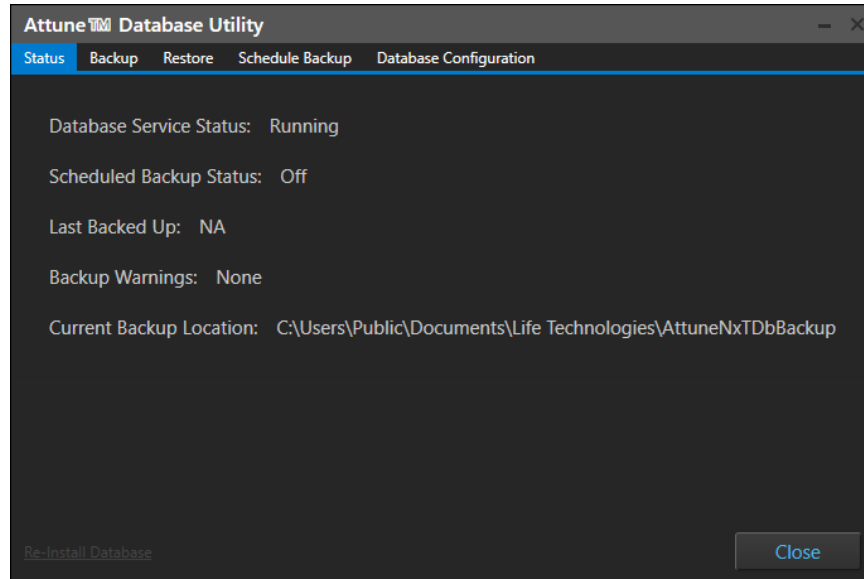
The visibility of the tabs is dependent on the user role. All users can see the **Status**, **Backup**, and **Restore** tabs. Only a **System Administrator** or **Administrator** can see the **Schedule Backup** and **Restore** tabs.

By default, **Status** tab is selected on logging into the application.



## Status tab

The **Status** tab displays the current status of the scheduled backup, the last backup time and date, the backup path, and any errors reported by the backup.



### Database Service Status

- The **Database Service Status** checks to see if the database service is running in the background. Typically, this displays the status of **Running**.
- If the backup service is **Stopped**, an option to start the backup service appears as the **Start Service** button next to the text. Click the **Start Service** button to start the backup.
- Successful start changes the status to **Running** and hides the **Start Service** button.

### Scheduled Backup Status

The **Scheduled Backup Status** shows the current **On** or **Off** state of the scheduled backup.

### Last Backed Up

- The **Last Backed Up** displays the date and time of the last backup in system locale format with the time shown in 24 hour format.
- If no backup has occurred, the status displays **NA**.

## Backup Warnings

The **Backup Warnings** displays any warnings or errors that have occurred when performing a scheduled backup.

If there are no warnings or errors, the status displays **None**. Possible errors include:

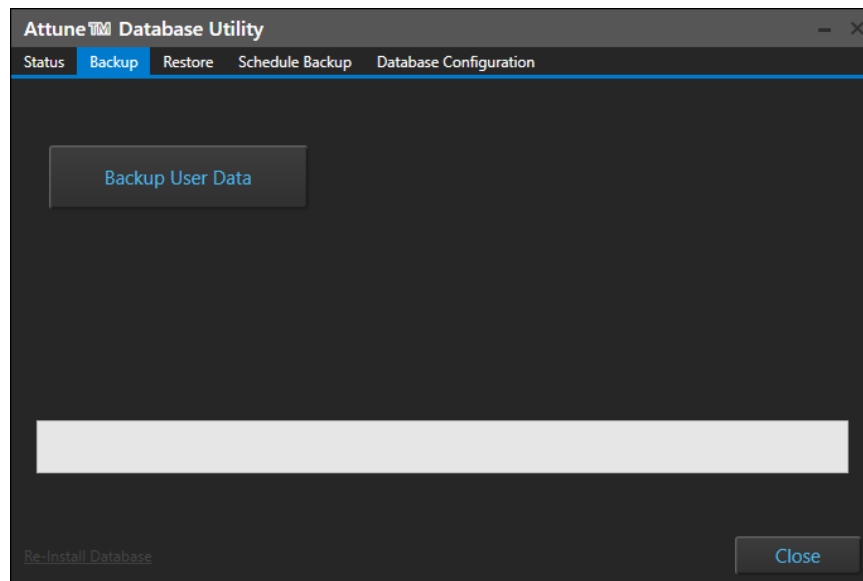
- **Insufficient Disk Space** warns that the attempted backup or database restore would consume > 90% of the free disk space.
- **Security Error** indicates that the user does not have permission to write files or folders at the destination.
- **Directory Not Found** indicates that the destination directories have been removed or renamed during backup or restore process.
- **Path Too Long** indicates that the file path name exceeds the number of characters allowed by the system.
- **Other exceptions** lists the last of the backup fails due to unforeseen problems.

## Current Backup Location

The **Current Backup Location** displays the current file path of the scheduled backup. If no path is specified, the status displays **NA**.

## Backup tab

The **Backup** tab enables the current user to back up all user data and the database to the location specified by the user (On demand backup of user data).



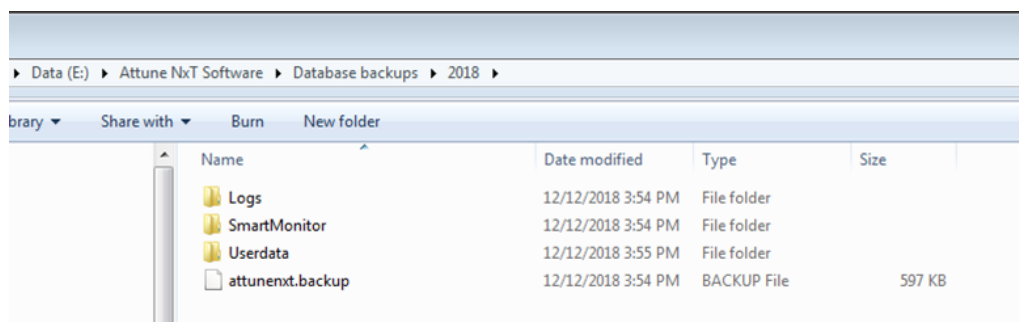


## Backup User Data (On demand backup of user data)

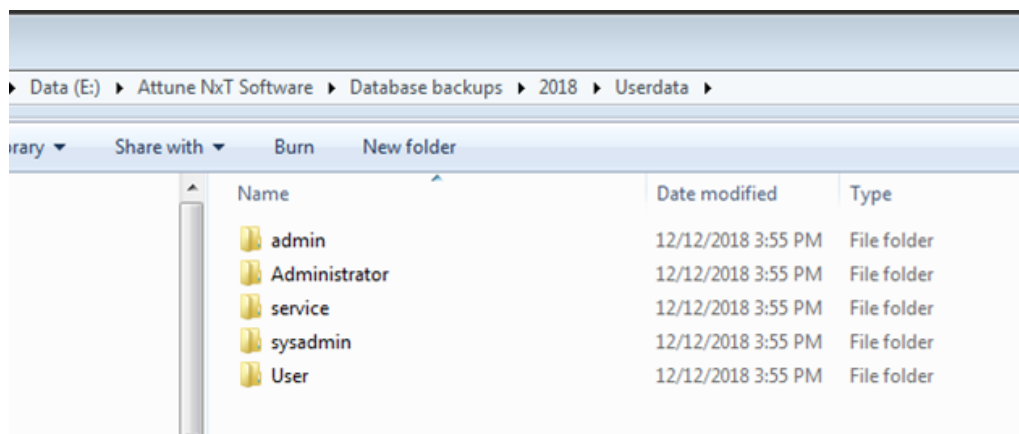
Click the **Backup User Data** button to open the file browser and select a location to back up the database and the corresponding data.

After confirming a location for the data backup, the database utility copies all user data in **\Users\Public\Public Documents\Thermo Fisher Scientific\AttuneNxt\Userdata**, then exports the database to the specified location.

- The database backup file that is created is called **attunenxt.backup** and includes information regarding the data structure.
- The data that are copied during the backup include copies of the Log files, Smart Monitor Data, and Userdata folders, which are in folders at the backup location.



- The Userdata folder contains subfolders for each account created in the software, including the default Service, System Administrator (sysadmin) and Administrator (admin) accounts, as well as custom, user-created accounts.



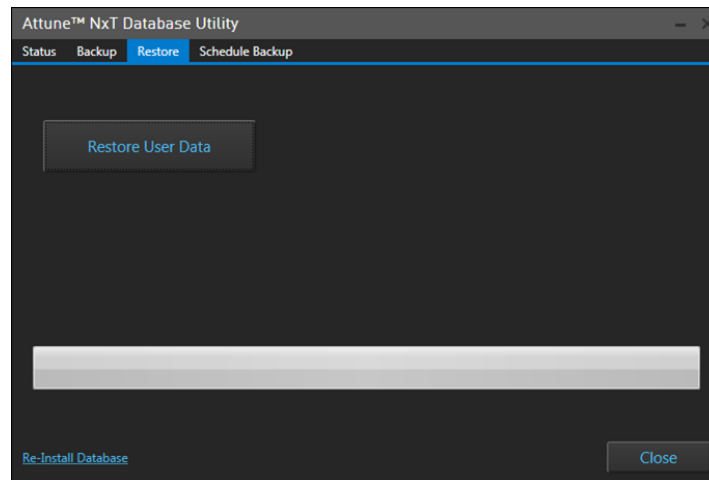
- In the folder for each user profile are data for each user account, including saved FCS files, and data associated with Experiment and Plate files, including Instrument settings, Workspaces, Compensation settings, Run protocols, Heat Map settings, Print settings, Overlay settings, Sample List settings, and Results Table settings.
- After a database backup has been completed once, if the same database backup location is selected for a secondary backup, only the files that have changed since the last backup are recopied. Any data files that have not changed are not recopied to the backup location. The new backup file replaces the previously saved **attunenxt.backup** file.

## Progress

A progress bar shows the progress of the backup operation as a percentage towards completion. The progress bar shows the current status as **Copying data** or **Backing up database**.

## Restore tab

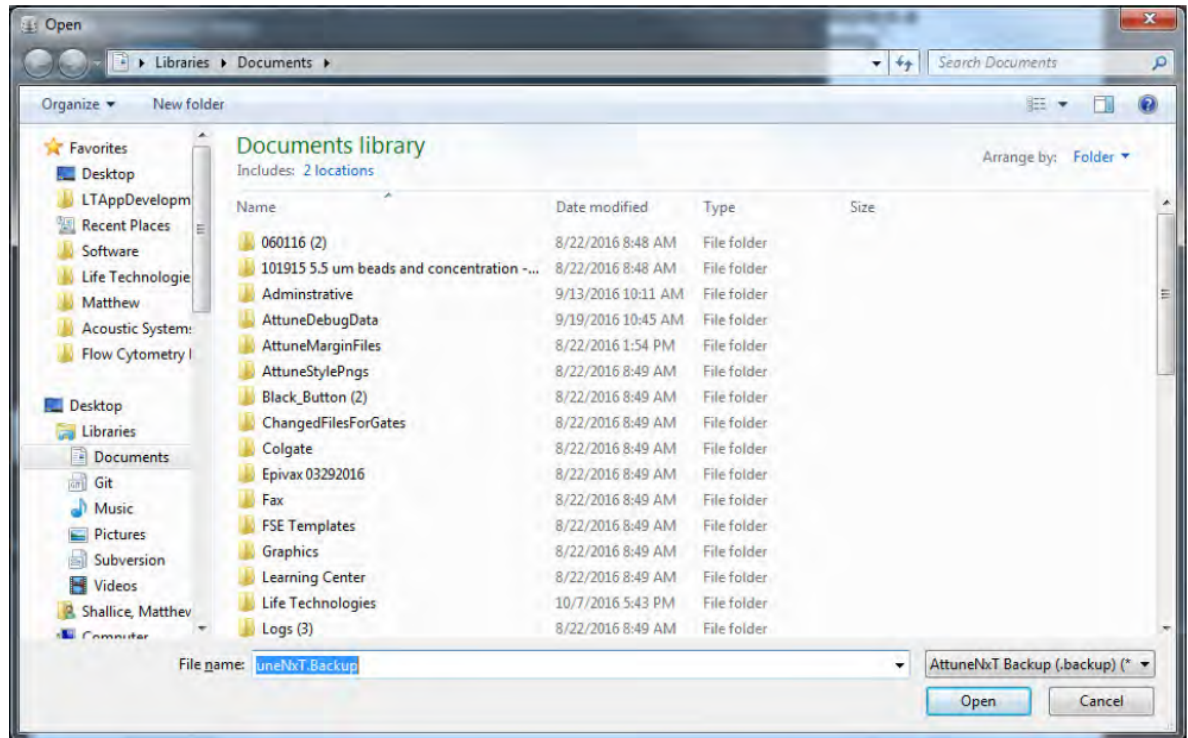
The **Restore** tab enables the user currently signed in to restore all user data and the database using a previously saved database.



## Restore User Data

**Restore User Data** operation cannot be completed when the Attune™ Cytometric Software is open. If the application is running, this option is disabled and displays a tooltip “This operation cannot be performed while the application is running.”

1. Click **Restore User Data** to open the **Open File dialog**:



2. Select a backup file and click **Open** to start the restore operation.

---

**IMPORTANT!** If the current database includes data that is not saved in the database backup file, new data are overwritten and lost as part of the database restoration process. To prevent data loss from occurring, export the data to a secondary location before initiating the database restoration.

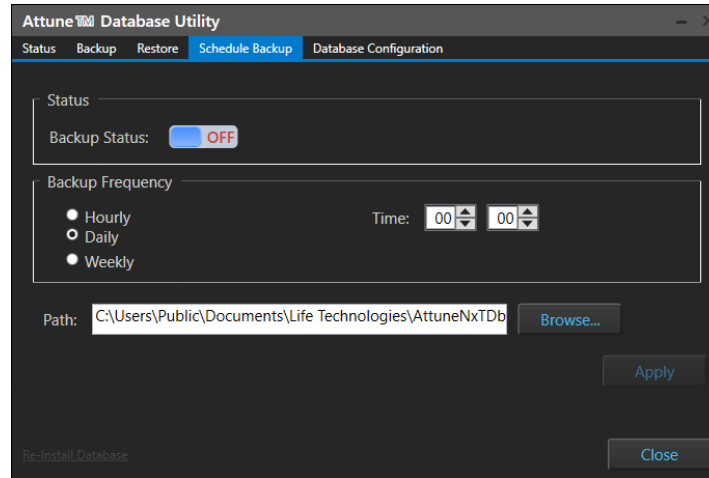
---

## Progress

A progress bar shows the progress of the restore operation as a percentage towards completion. The progress bar shows the current status as **Copying data** or **Restoring database**.

## Schedule Backup tab

The **Schedule Backup** tab enables an authorized user to set the backup policy and schedule. This tab is only visible to the **System Administrator** or **Administrators**.



### Status

The **Status** group includes a toggle control to turn the backup on or off.

- By default the **Backup Status** option is set to **OFF**.
- When turned **ON**, the utility attempts to backup to the specified path and frequency.
- If an error occurs when performing a scheduled backup, the error is displayed on the **Status** tab (“Status tab” on page 817).
- When turned **OFF**, no automatic backup occurs.
- If a valid path is not entered, a warning message is displayed when the user tries to turn the status **ON**: “Valid path is required to schedule an automatic backup.”

### Backup Frequency

The **Backup Frequency** sets the frequency of the scheduled backup. You can set the scheduled backup to run **Hourly**, **Daily**, or **Weekly**.

- By default, the **Backup Frequency** is set to **Daily**.
- **Time** specifies the time to start the backup. The time format is in 24 hour time as HH:MM.
- By default, the backup is scheduled to run at **00:00**.
- When the **Hourly** option is selected, the **Time** control is not visible and the backup runs at the start of the hour (:00).
- When the **Weekly** option is selected, the **Day** must be specified. By default, the day is set to **Sunday**.

## Path

**Path** specifies the file path of the backup folder. You can type the file path of the backup folder directly into the **Path** textbox or click **Browse** to select it from the **File path dialog**.

If the file path does not exist, the folder is automatically created when the application is closed.

## Copied files

The first time the database backup is completed, all data are copied to the backup location and an **attunenxt.backup** file is created.

Every time the scheduled backup occurs thereafter, only data that has changed after the last backup are copied; any data files that have not changed are not recopied to the backup location.

The new backup file replaces the previously saved **attunenxt.backup** file each time the scheduled backup is completed.

## Reinstall Database

The **Reinstall Database** hyperlink is only visible to the **System Administrator** and **Administrators**. Reinstallation of the database replaces the existing database with a new, blank database.

---

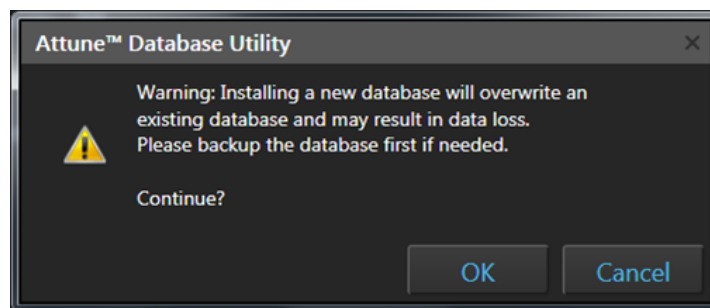
**Note:** Database reinstallation completely erases all data in the current database. However, it does not erase data saved in the previous backup.

---

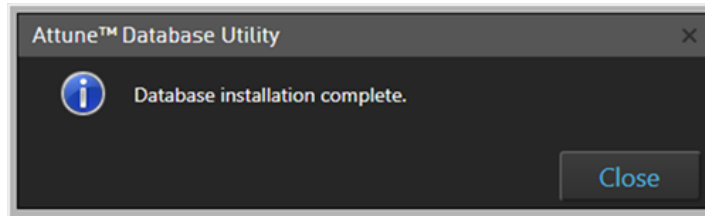
**IMPORTANT!** If the backup scheduler is set to back up user data at a specified time, the first scheduled backup that occurs after a database reinstallation overwrites all previously saved data. To prevent data loss from occurring, copy the **Userdata folders** and **attunenxt.backup** file to an alternate location before using the scheduled backup procedure.

---

1. Click the **Reinstall Database** hyperlink. A confirmation dialog opens.



- Click **OK** to reinstall the Attune™ Database. When the reinstallation is completed, the following confirmation dialog appears:




---

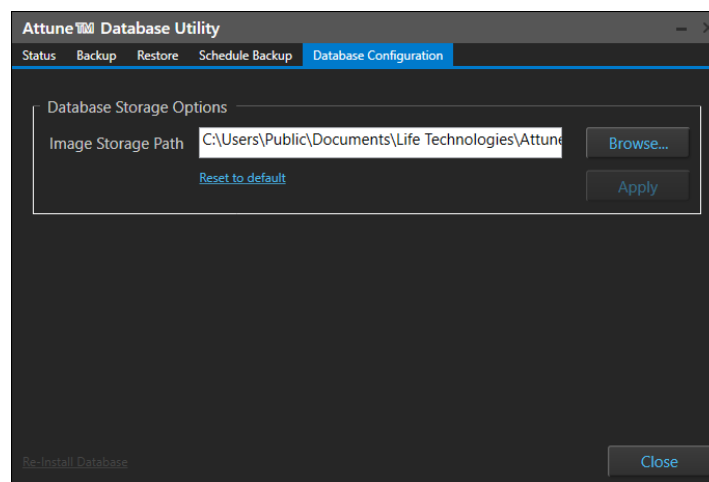
**Note:** If the Attune™ Cytometric Software is running, the **Reinstall Database** hyperlink is disabled and displays a tooltip “This operation cannot be performed while the application is running.”

---

## Database Configuration tab

The **Database Configuration** tab enables an authorized user to set the **Database Storage Options** by selecting the **Image Storage Path** for images captured with the Attune™ CytPix™ Flow Cytometer.

This tab is only visible to the **System Administrator** or **Administrators** when an Attune™ CytPix™ Flow Cytometer is in use.



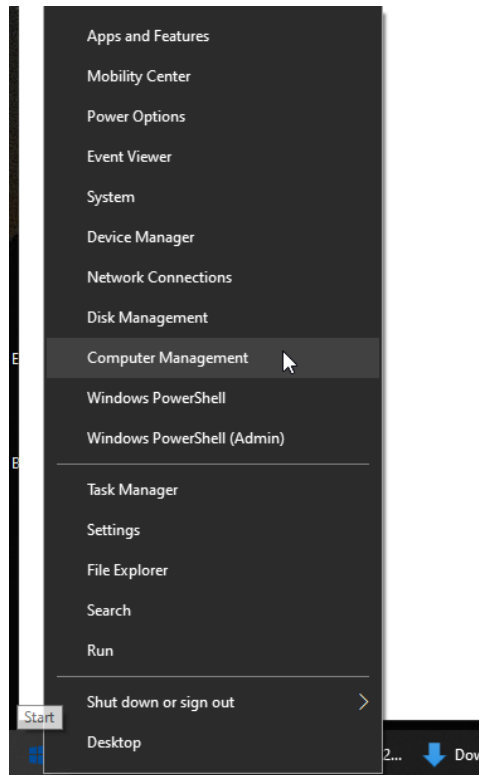
## Image Storage Path

- You can type the File Path of the backup folder directly into the **Image Storage Path** textbox or click **Browse** to select it from the **File path dialog**.
- If the File Path does not exist, the folder is automatically created when the application is closed.
- The default image storage path is:  
**C:\Users\Public\Documents\LifeTechnologies\AttuneCytPix\_Images**
- To reset the image storage path, click **Reset to default**.

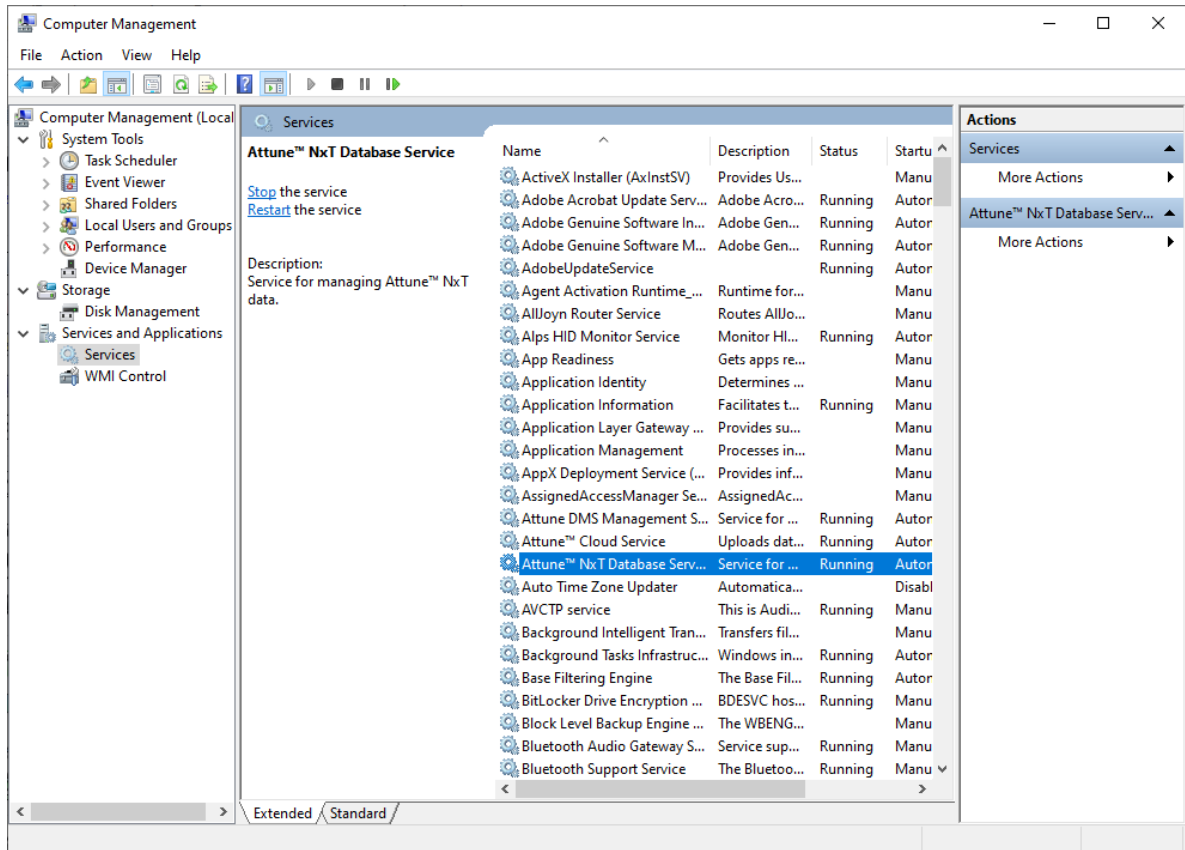
## Attune™ Database Service

The **Attune™ Database Service** is a Windows™ data backup service that runs in the background and performs the backup at the scheduled time. You can manage the service via the **Windows™ Management Console**.

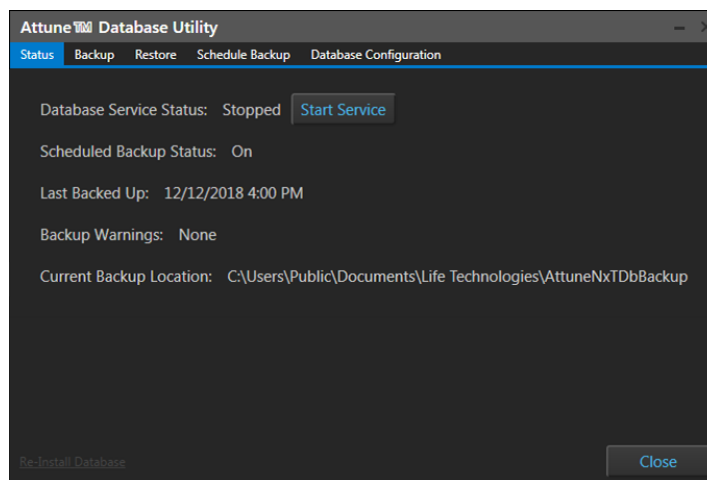
1. To access the **Windows™ Management Console**, right-click the **Start** icon on the computer desktop, then select **Computer Management**.



- In the **Computer Management** dialog, under **Services and Application**, select **Services**, then select **Attune™ NxT Database Service**.



- You can start or stop the service from the **Windows™ Management Console** or double-click the service to get more options.
- After the service is stopped, you can start the service again from the **Attune™ Database Utility** application.







# FCS file reference

## Overview

This chapter describes the FCS keywords that are used when generating FCS files from data acquired from the Attune™ NxT Flow Cytometer.

- When in the Record mode, the FCS is recorded regardless of whether the stop criteria are met.
- For appended data sets, the time and event offsets (elapsed time from \$BTIM in ms and \$TOT of previous dataset; see below) are added to the time and event parameter values (when enabled) in the new data stream.
- The FCS files generated by the Attune™ Cytometric Software when recording or exporting conform to the FCS standard as described by the Data File Standards Committee of the International Society for Analytical Cytology (ISAC) for FCS 3.1.

## FCS HEADER format

The *FCS HEADER* conforms to the FCS standard as described by the Data File Standards Committee of the International Society for Analytical Cytology (ISAC). An example table is displayed below.

Contents	Start by position	End byte position
FCS 3.1	00	05
ASCII(32) – space characters	06	09
ASCII-encoded offset to first byte of TEXT segment	10	17
ASCII-encoded offset to last byte of TEXT segment	18	25
ASCII-encoded offset to first byte of DATA segment	26	33
ASCII-encoded offset to last byte of DATA segment	34	41
ASCII-encoded offset to first byte of ANALYSIS segment	42	49
ASCII-encoded offset to last byte of ANALYSIS segment	50	57
ASCII-encoded offset to user-defined OTHER segment	58	beginning of next segment

## FCS TEXT segment

- The delimiter used in the TEXT segment of the FCS file is / (ASCII 47). If the delimiter appears in a keyword or keyword value, it is immediately followed by a second delimiter //.
- For custom keywords, the # symbol precedes the keyword name instead of the \$ symbol.
- The FCS Keywords in the FCS TEXT segment are listed in alphabetical order.
- The parameter description keywords (\$PnN, \$PnR, etc.) in the FCS TEXT segment are numbered consecutively and are listed in the order the parameters are written to the file.

### Required keywords

Keyword	Description	Source of Value
\$BEGINANALYSIS	Byte-offset to the beginning of the ANALYSIS segment	This is set to "0"
\$BEGINDATA	Byte-offset to the beginning of the DATA segment	By default, this is set at 8192. If the TEXT segment exceeds this byte offset, the value will be set to ensure correct alignment of the DATA segment. This byte-offset is a duplicate of the offset contained in the HEADER. If the DATA segment exceeds the first 99,999,999 bytes, the offset in the HEADER will be set to 0 and this value will take precedence.
\$BEGINSTEXT	Byte-offset to the beginning of a supplemental TEXT segment	This is set to "0"
\$BYTEORD	Byte order for data acquisition computer	This is set to "1,2,3,4"
\$DATATYPE	Type of data in DATA segment (ASCII, integer, floating point)	This is set to "F"
\$ENDANALYSIS	Byte-offset to the end of the ANALYSIS segment	This is set to "0"
\$ENDDATA	Byte-offset to the end of the DATA segment	This is dynamically updated based on the offset of the last DATA byte.
\$ENDSTEXT	Byte-offset to the end of a supplemental TEXT segment	This is set to "0"
\$MODE	Data mode (list mode, histogram)	This is set to "L"
\$NEXTDATA	Byte offset to next data set in the file	This is set to "0"
\$PAR	Number of parameters in an event	From the total number of enabled parameters, including custom parameters, recorded in to the FCS file.
\$PnB	Number of bits reserved for parameter number n	This is set to "32"

\$PnE	Amplification type for parameter n	This is set to 0,0
\$PnR	Range for parameter number n	By default this is set to $2^{26}$ for Event. <sup>[1]</sup> By default this is set to $2^{26}$ for Time.* By default this is set to $2^{20}$ for all Height measurements.* This is set to $2^{20}$ for all Area measurements.* By default this is set to $2^{10}$ for all Width measurements.*
\$TOT	Total number of events in the data set	Number of events in the file (updated when appended).

<sup>[1]</sup> Data values transferred from the instrument are constrained to be within the \$PnR limits.

## Optional keywords

Keyword	Description	Source of Value
\$ABRT	Events lost due to data acquisition electronic coincidence	From datastream when “Exclude coincident events” is selected in the Instrument Settings Panel as described on “Exclude coincident events check box” on page 395.
\$BTIM	Clock time at beginning of data acquisition	The system clock time at the time the recorded data stream commences.
\$CELLS	Description of objects measured	From Sample information Notes field as described on “Sample information” on page 473.
\$COM	Comment	
\$CYT	Type of flow cytometer	From the Attune™ model name and number as described on “About Attune™ cytometric software” on page 702.
\$CYTSN	Flow cytometer serial number	From the cytometer serial number as described on “About Attune™ cytometric software” on page 702.
\$DATE	Date of data set acquisition	The system clock date at the time the recorded data stream commences.
\$ETIM	Clock time at end of data acquisition	The system clock time at the time the recorded data stream ends. For appended data, this is based on the time of the last addition to the data file.
\$EXP	Name of investigator initiating the experiment	From the name of the PI, if specified, for the currently logged in user as specified on “Add or edit an account” on page 705.
\$FIL	Name of the data file containing the data set	From the sample name at the time the data is recorded. This is the pathless filename and extension.
\$INST	Institution at which data was acquired	From the institution name field as specified on “Add or edit an account” on page 705.

(continued)

Keyword	Description	Source of Value
\$LAST_MODIFIED	Timestamp of the last modification of the data set	The time a data file is appended or updated based on exporting the dataset.  This is only added when the \$ORIGINALITY is not set to Original.
\$LAST_MODIFIER	Name of the person performing last modification of a data set	The name of the currently logged in user when a data file is appended or updated based on exporting the dataset.  This is only added when the \$ORIGINALITY is not set to Original.
\$LOST	Number of events lost due to computer busy	From the data stream
\$OP	Name of flow cytometry operator	From the name (first and last) of the currently logged in user
\$ORIGINALITY	Information whether the FCS data set has been modified (any part of it)	For appended datasets, the keyword value is set to "Appended".  For exported data where the TEXT segment or data segment is modified, the keyword value is set to "DataModified".  When the dataset has not been modified, this is set to "Original".
\$PLATEID	Plate identifier	From the "Plate ID" field s. If it is left blank, the keyword is the plate name.
\$PLATENAME	Plate name	From the "Plate Name" field.
\$PnF	Name of optical filter for parameter n	From the instrument configuration information for the specified detector.
\$PnL	Excitation wavelength(s) for parameter n	From the instrument configuration information for the specified detector.
\$PnS	Name used for parameter n	From the concatenation of the "target and label". When absent, the \$PnS is equivalent to the \$PnN.
\$PnV	Detector voltage for parameter n	From the voltage for each detector at the time the file was recorded.
\$PnN	Channel name for parameter n	From the channel name for the specified detector.
\$PROJ	Name of the experiment project	From the experiment name.
\$SMNO	Specimen (e.g., tube) label	From the group name.

(continued)

Keyword	Description	Source of Value
\$SPILLOVER	Fluorescence spillover matrix	From the Experiment-level spillover matrix state at the time the file is recorded or exported.
\$SRC	Source of the sample (sample name, cell types)	
\$SYS	Type of computer and its operating system	From the system attributes including computer type and operating system.
\$TIMESTEP	Time step for time parameter	This will be set to 0.001 but should be provided by the instrument API.
\$TR	Trigger parameter and its threshold	From the threshold settings set at the time a file was recorded in the format (Boolean_Channel, Boolean_Channel...).
\$VOL	Volume of sample run during data acquisition	From the instrument API. Volume is updated to reflect total volume for all draws of a multidraw, and, if appended, the volume is totaled.
\$WELLID	Well identifier	From the well location (i.e., A1, A2,...) for a Plate sample.  For tubes recorded in a Plate Experiment, the tube location for the Tube sample is recorded (i.e., T1, T2...).

## Custom keywords

Keyword	Description	Source of Value
#TOTALVOLUME	Volume of sample run during data acquisition	From the instrument API. Volume is updated to reflect total volume for all draws of a multidraw, and, if appended, the volume is totaled.
#PnTarget	The target name for parameter n	From the target field for each parameter within the Instrument Settings Panel.
#PnLabel	The label name for parameter n	From the label field for each parameter within the Instrument Settings Panel.
#FLOWRATE	The flow rate of the sample during data acquisition	From the instrument API or collection panel state.
#LASER1ASF	Area Scaling Factor used for channels recorded off of laser 1	Area scaling factor for detectors off of laser 1 at time of recording.
#LASER2ASF	Area Scaling Factor used for channels recorded off of laser 2	Area scaling factor for detectors off of laser 2 at time of recording.

(continued)

Keyword	Description	Source of Value
#LASER3ASF	Area Scaling Factor used for channels recorded off of laser 3	Area scaling factor for detectors off of laser 3 at time of recording.
#LASER4ASF	Area Scaling Factor used for channels recorded off of laser 4	Area scaling factor for detectors off of laser 4 at time of recording.
#LASER1DELAY	Laser Delay of Laser 1	The laser delay setting for laser 1 at time of recording. Only recorded if laser 1 is present and turned on.
#LASER2DELAY	Laser Delay of Laser 2	The laser delay setting for laser 2 at time of recording. Only recorded if laser 2 is present and turned on.
#LASER3DELAY	Laser Delay of Laser 3	The laser delay setting for laser 3 at time of recording. Only recorded if laser 3 is present and turned on.
#LASER4DELAY	Laser Delay of Laser 4	The laser delay setting for laser 4 at time of recording. Only recorded if laser 4 is present and turned on.
#WINEXT	Window Extension	The window extension setting at time file is recorded
#PTRESULT	Performance Test Result	Performance Test status at time of recording (Pass/Fail)
#PTDATE	Date of most recent performance test	Date of most recent performance test at time of recording.
#WIDTH THRESHOLD	Width Threshold	The width threshold setting at the time of recording. Only needed, if width is a recorded measurement.
#LASER1COLOR	Laser 1 Color	Laser 1 Color at time of recording. Specified in terms of wavelength in units of $\mu\text{m}$ .
#LASER2COLOR	Laser 2 Color	Laser 2 Color at time of recording. Specified in terms of wavelength in units of $\mu\text{m}$ .
#LASER3COLOR	Laser 3 Color	Laser 3 Color at time of recording. Specified in terms of wavelength in units of $\mu\text{m}$ .
#LASER4COLOR	Laser 4 Color	Laser 4 Color at time of recording. Specified in terms of wavelength in units of $\mu\text{m}$ .
#LASERCONFIG	The laser configuration of the Attune™ Instrument	Laser configuration at time of file recording.

(continued)

Keyword	Description	Source of Value
#DRAWnVOLUME	Acquisition volume of the n <sup>th</sup> draw in a multi-draw	Instrument API will provide this value upon completion of nth draw. Only recorded if the TIME acquisition parameter is enabled.
#DRAWnSTART	Acquisition start time in terms of the TIME parameter	Instrument API will provide this value upon start of nth draw. Only recorded if the TIME acquisition parameter is enabled.
#DRAWnSTOP	Acquisition stop time in terms of the TIME parameter	Instrument API will provide this value upon completion of nth draw. Only recorded if the TIME acquisition parameter is enabled.

## Storage of user-defined keywords in the FCS file

User-defined keywords are stored as custom keywords when recording FCS files.

- When recording FCS files, the keyword values are written to the FCS file as customer keyword value pair, where the # symbol is added to the front of the name to indicate that it is a custom keyword.
- The custom keywords are also considered when exporting an FCS file and the user is provided with the option to update these keywords.

## Storage of compensation values in the FCS file

- If an Experiment Compensation exists prior to recording a sample, the spillover values are recorded into the FCS file using the \$Spillover keyword.
- Only the parameters that are included in the compensation calculation are included in the \$Spillover keyword value.
- Compensation controls only contain the identity matrix.



# Attune™ Cytometric Software image processing parameters

## Image processing parameters

Feature	Description	PnR <sup>[1]</sup>
<b>Intensity and Texture features</b>		
AverageIntensity	Average intensity of all pixels within an object	2 <sup>10</sup>
AverageNormIntensity	$100 \times \text{AverageIntensity} / (0.5 \times 2^{(\text{BitsPerPixel} - 1)})$	
CVIntensity	$100 \times \text{StandardDeviationIntensity} / \text{AverageIntensity}$	
CVNormIntensity	$100 \times \text{StandardDeviationNormIntensity} / \text{AverageNormIntensity}$	
EntropyIntensity	Entropy of intensity distribution of all pixels within an object	
KurtosisIntensity	Kurtosis of intensity distribution of all pixels within an object <sup>[2]</sup>	
MaxIntensity	Maximum intensity of all pixels within an object	2 <sup>10</sup>
MinIntensity	Minimum intensity of all pixels within an object	2 <sup>10</sup>
SkewnessIntensity	Skewness of intensity distribution of all pixels within an object <sup>[3]</sup>	
StandardDeviationIntensity	Standard deviation of intensity of all pixels within an object	2 <sup>10</sup>
StandardDeviationNormIntensity	$100 \times \text{StandardDeviationIntensity} / (0.5 \times 2^{(\text{BitsPerPixel} - 1)})$	
TotalIntensity	Total intensity of all pixels within an object	
<b>Object features</b>		
ParticleCount	Number of cells within the identified object	2 <sup>3</sup>
<b>Pixel features</b>		
NumPixels	Number of pixels contained within identified objects	2 <sup>16</sup>
<b>Shape features</b>		
AreaSquareMicrons	Area of the object measured within mask based on pixel count = NumPixels × (PixelSize.MicronsX × PixelSize.MicronsY)	2 <sup>13</sup>
CircularityPercent	Percent circularity of an object = 100 / PerimeterToArea (100 for a circular object)	2 <sup>9</sup>



(continued)

Feature	Description	PnR <sup>[1]</sup>
EccentricityPercent	Eccentricity of an ellipse = $100 \times \text{Sqrt}(1 - \text{ShortAxisMicrons}^2 / \text{LongAxisMicrons}^2)$	0 to 100
MajorDiameterMicrons	Distance across an ellipse along its long axis = $\text{MajorRadius} \times 2 \times \text{PixelSize.MicronsX}$	2 <sup>7</sup>
MinorDiameterMicrons	Distance across an ellipse along its short axis = $\text{MajorRadius} \times 2 \times \text{PixelSize.MicronsX}$	2 <sup>7</sup>
PerimeterMicrons	Perimeter of an object	2 <sup>15</sup>
PseudoDiameterMicrons	Diameter of a circle with an area equal to the area of the object = $2 \times \text{Sqrt}(\text{Object.Area in } \mu\text{m}^2 / \text{pi})$	2 <sup>15</sup>
MinorMajorRatioPercent	Short to long axes ratio of an object as a percentage = $100 \times \text{ShortAxisMicrons} / \text{LongAxisMicrons}$	0 to 100
<b>System features</b>		
ConfidenceScore	Indicates that one or more objects intersects with the FOV of the image	
IsOnBorder	Indicates that one or more objects intersects with the FOV of the image	
IsProcessable	Indicates that the image is processable	
IsProcessed	Indicates that the image was processed. This is generated when the image processing FCS file is loaded and merged with the raw FCS file data.	

[1] \$PnR is the range for the selected parameter n.

[2] Kurtosis of intensity measures the peakedness of the distribution of all pixels within an object. See “Kurtosis” on page 836.

[3] Skewness of intensity measures the degree of asymmetry in the pixel data of an object. See “Skewness” on page 836.

## Kurtosis

Kurtosis is measures how peaked a histogram is and it is based on the size of a distribution's tails. The kurtosis of a normal distribution is 0. Distributions with short tails compared to a normal distribution have negative kurtosis (platykurtic) and distributions with relatively long tails have positive kurtosis (leptokurtic).

In the context of image processing, kurtosis describes whether distribution of gray tones is more spread-out (flat) or more concentrated around the mean (peaked).

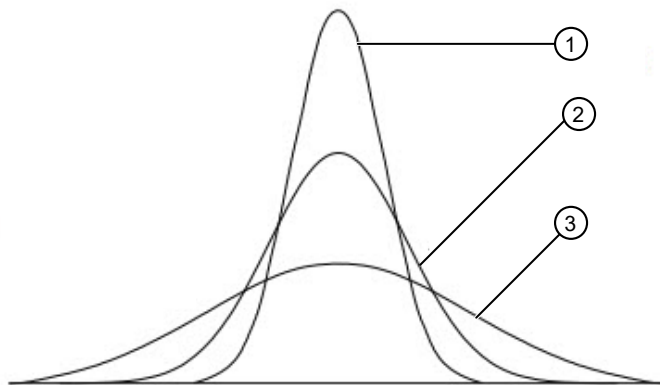


Figure 140 Kurtosis

- ① Positive kurtosis (leptokurtic)
- ② Normal distribution (mesokurtic)
- ③ Negative kurtosis (platykurtic)

## Skewness

Skewness measures the degree of asymmetry exhibited by the data. If skewness equals zero, the histogram is symmetric about the mean.

In the context of image processing, skewness indicates the imbalance between the number of pixels that are darker or brighter than the mean.

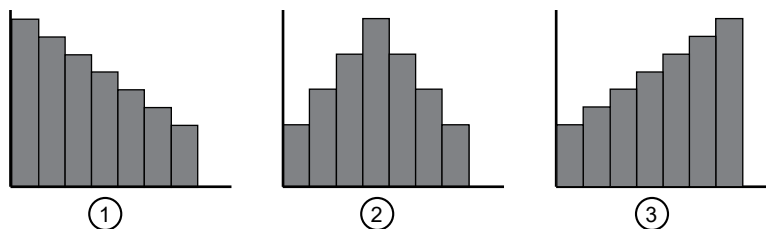


Figure 141 Skewness

- ① Positively skewed histogram
- ② Symmetric distribution histogram
- ③ Negatively skewed histogram



# Data management in Attune™ Cytometric Software

## Overview

This section describes options for managing (saving and/or deleting) data in the Attune™ Cytometric Software. Note that it is critical for you to review all instructions completely before executing the steps outlined in this section. We recommend that you archive data before deletion.

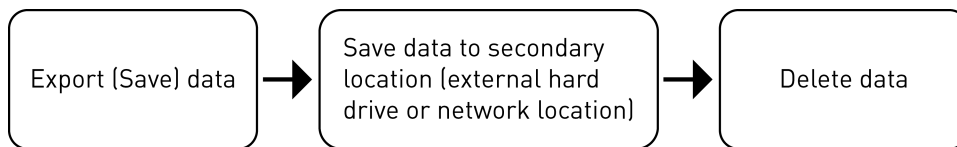


Figure 142 Data management workflow

## Contents

- Saving user data
- Types of data included in the Attune™ Cytometric Software
- Disk usage
- Data storage warnings
- Change low disk space warning options
- Choosing one or more data archiving methods
- Database backup
- How to save and export Plate and Tube Experiment files
- How to save and export FCS files
- How to save and export FCS and image data
- How to save and export images
- How to use the Attune™ Database Utility program
- How to restore data using a saved database
- How to copy a database to a different computer
- Deletion of Experiment and Plate files
- Reinstallation of a new database
- Deleting a user account to remove data
- Permanent removal of old Baseline and Performance Test data



## Types of data in Attune™ Cytometric Software

There are many types of data associated with the Attune™ instrument and software (Figure 143). Users create “data” each time they create a Tube or Plate Experiment and when FCS or image files are recorded during sample acquisition. Data also refers to other inputs into the software, such as instrument management information that allow the cytometer to collect samples, retain performance data and log information. User management data are retained by the software and provides instructions to manage user accounts, security, and user preferences. Data saved from analysis of results include information stored in the Heat Map, Results table, Overlay settings, and Image View, as well as image processing data.

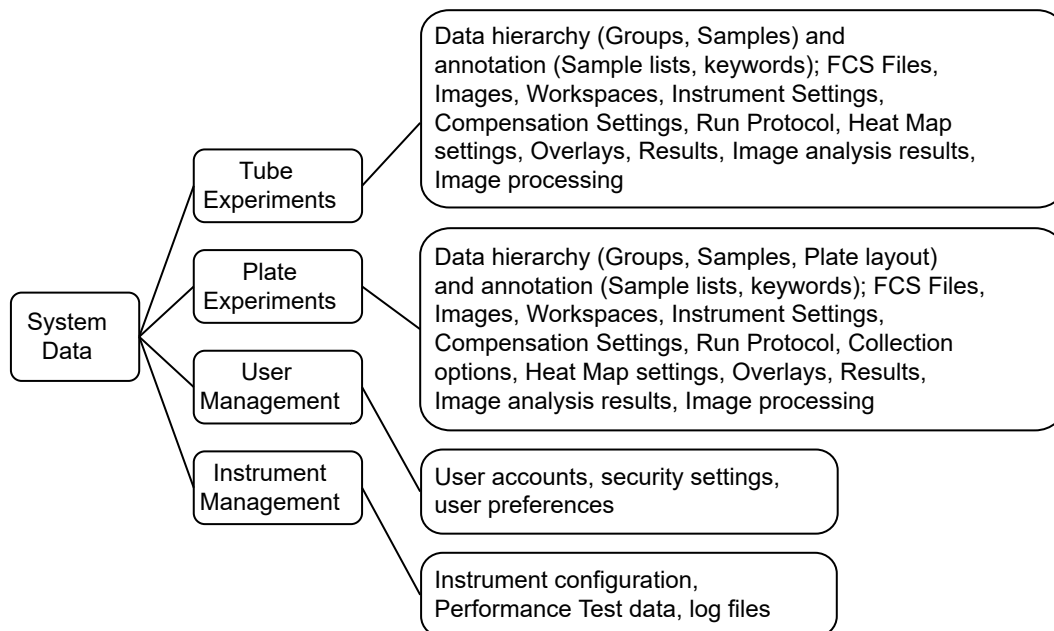
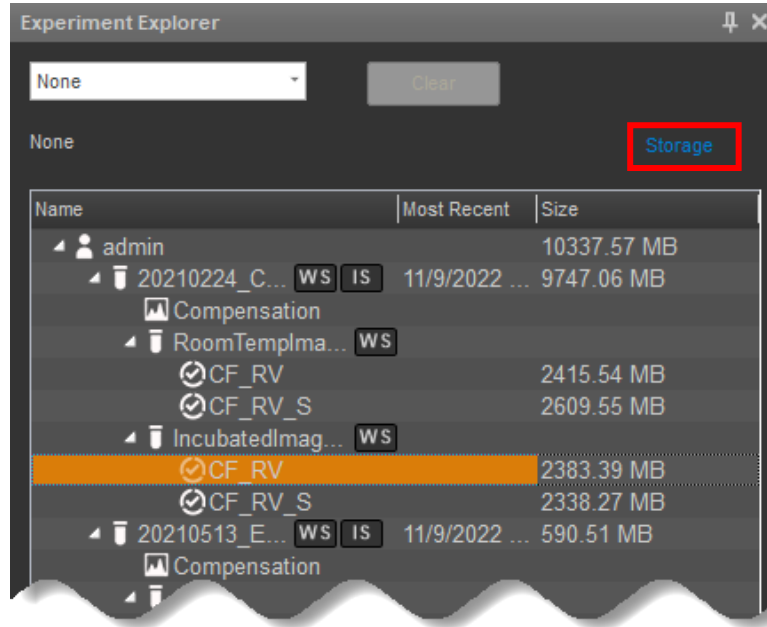


Figure 143 Data in the Attune™ Cytometric Software



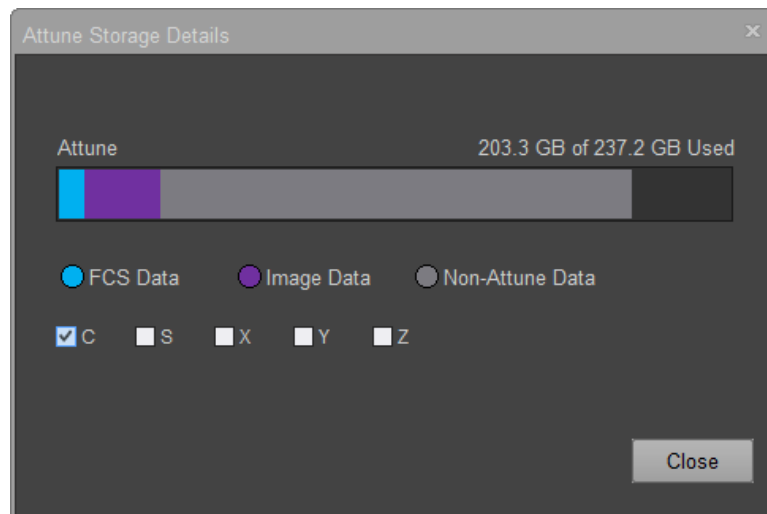
## Disk usage

To view the data storage details on the instrument, click the **Storage** hyperlink at the top of the **Experiment Explorer**, which opens the **Attune Storage Details** window.



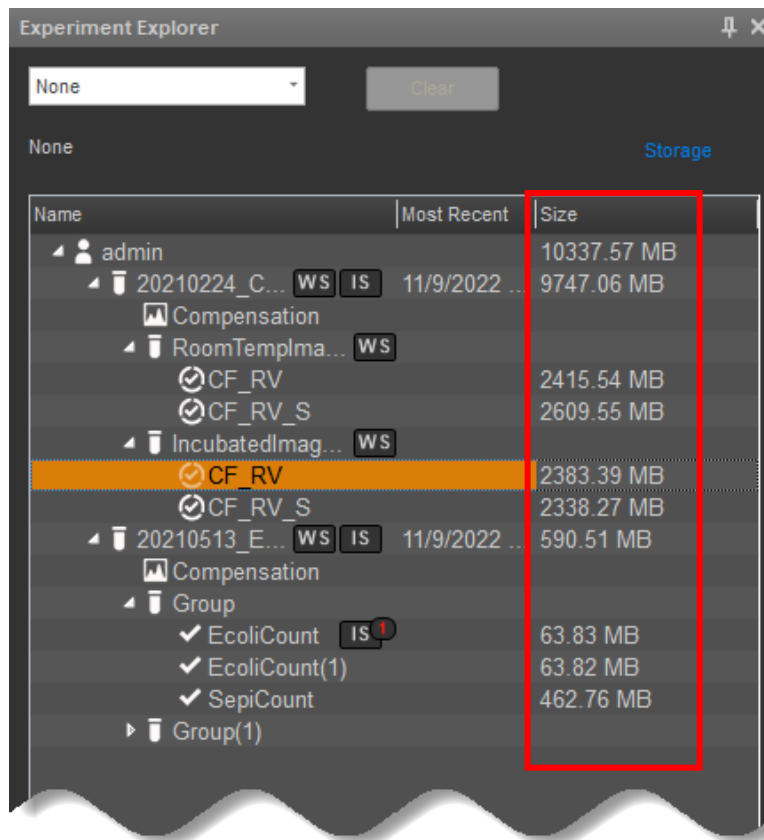
The **Attune Storage Details** displays the total used space and the total combined space broken into:

- FCS Data
- Image Data
- Non-Attune Data



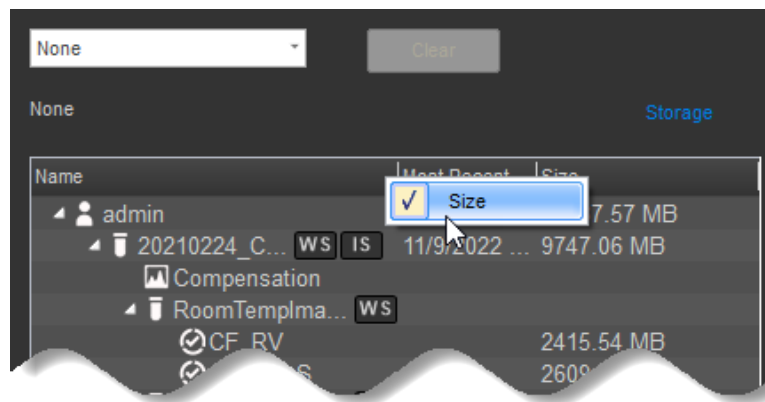


The size details for Experiments, Groups, and Samples are shown in the **Size** column in **Experiment Explorer**, which are calculated on sign in and when any changes are made to an experiment (data added or removed).



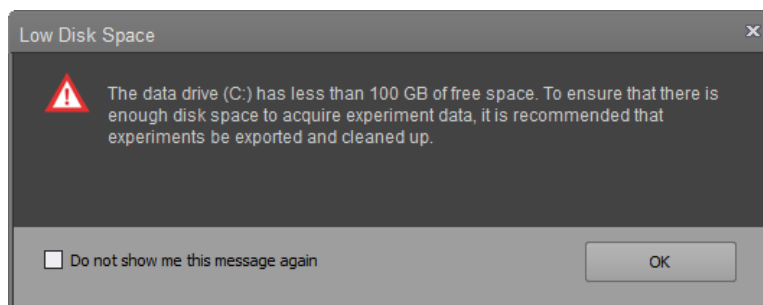


- The size data are displayed for all samples that contain data and include the size of the FCS file, the image data, and all metadata for the sample.
- The size data are displayed for the Experiment and include the summed size of all Sample data and any Experiment metadata.
- To sort the size data in ascending or descending order, click the **Size column header**.
- To hide or show the size column, right-click the **Experiment Explorer heading row**, then deselect or select **Size**.



## Data storage warnings

By default, when activating experiments, the Attune™ Cytometric Software displays the **Low Disk Space** warning if the available disk space is less than 100 GB on the primary data (image data and experiment data) drives.



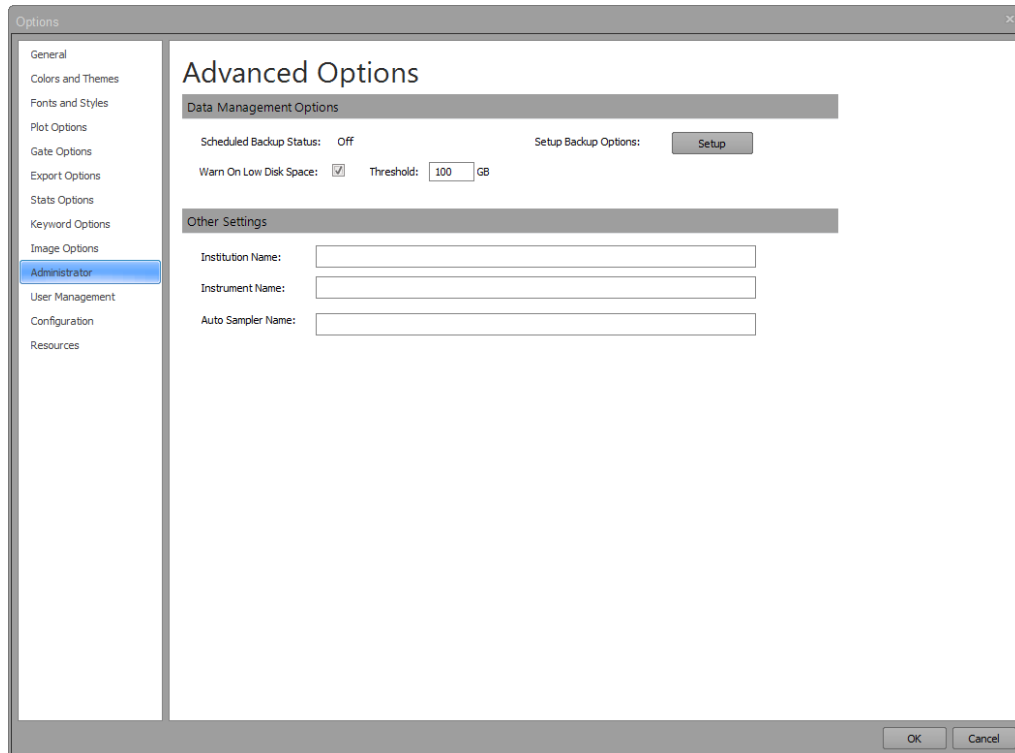
- To ignore and suppress the warning for the length of the session, select **Do not show me this message again**, then click **OK**.
- To disable the warning or to change the low disk threshold, see “Change low disk space warning options” on page 842.



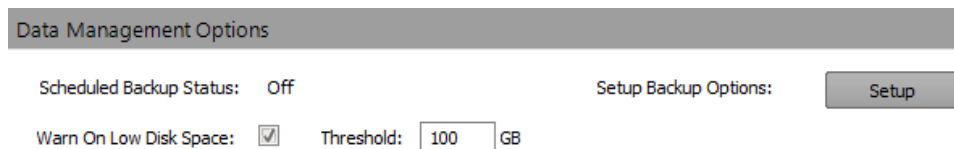
## Change low disk space warning options

Users with Administrator or System Administrator accounts can change backup and low disk space warning options in the **Administrator** tab of the **Options** dialog.

1. On the **Quick Access** toolbar, click **Options** to open the Options dialog, then select the **Administrator** tab.



2. To disable the Low Disk Space warning, deselect **Warn On Low Disk Space**.



3. To change the threshold of available disk space on the primary disk drives below which the Low Disk Space warning is displayed, enter the desired value (in GB) in the **Threshold** field.

---

**Note:** By default, the threshold is set to 100 GB. You can enter a threshold value of between 1 GB and 1000 GB.

---





## Options for saving user data

There are many options for saving user data from the Attune™ Cytometric Software. You can export Tube and Plate experiment files (.apx and .atx), individual FCS files (.fcs), FCS and image processing data, and image files from the software and store them in a secondary location. Alternatively, you can use the Attune™ Database Utility program to create a copy of all the experiments and associated data files in the software at the time the backup is created.

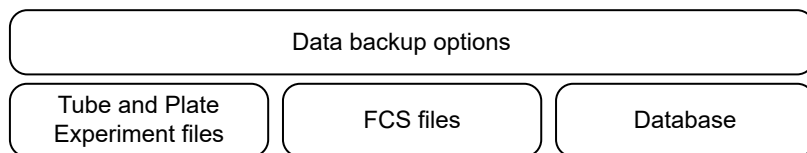


Figure 144 Options of saving user data from the Attune™ Cytometric Software

## Choosing one or more data archiving methods

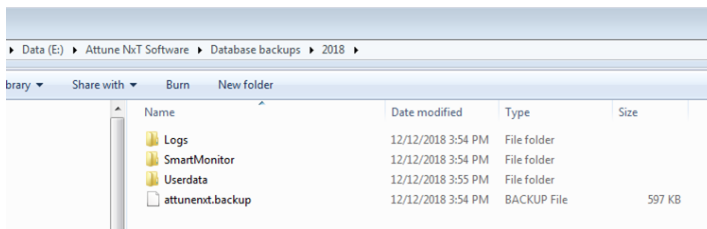
- When you export a Tube Experiment file as an .atx file, the information saved includes all data associated with the Experiment: Groups and Samples, recorded FCS files, Run Protocol settings, Instrument settings, Workspace settings, Compensation files and settings, Heat Map settings, Sample lists, Overlays, Results table settings, and print settings.
- When you export a Plate Experiment as an .apx file, the information that is saved includes the same information as an Experiment file, but also includes Plate layout information and the Plate collection settings.
- When you export FCS files, you can save the file as an FCS 3.0 or FCS 3.1 file data standard. The information that is saved includes the raw data acquired for the sample and associated sample information that is annotated using FCS file keywords (see “Appendix A, “FCS file reference”). Some keywords provide annotation and standardization about the sample (such as cytometer used for acquisition, time of acquisition, number of events, etc.), while others are Attune™ Cytometric Software custom keywords (including flow rate, etc.) or user defined custom keywords (such as dilution factor, sample ID, etc.). You can reimport FCS files individually into the Attune™ Cytometric Software or you can analyze them using other flow cytometry analysis software.
- When the Attune™ Database Utility program is used as a file backup system (see “Database backup” on page 844), the information that is saved includes all data files used in the software as described above and also includes user management settings (profiles, user settings, security settings), instrument configuration settings, baseline and performance test data, and instrument data logs.



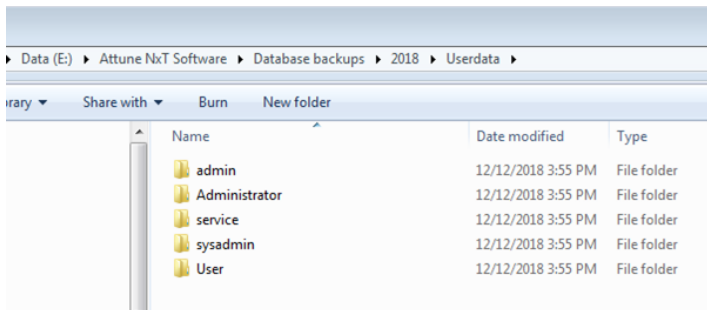
## Database backup

The Attune™ Database Utility (Chapter 29, “Attune™ Database Utility”) is a data backup program that is provided with the Attune™ Cytometric Software. This utility program enables you to back up user data to a custom location, either on-demand or at a specified, scheduled frequency.

- The database utility copies all user data in the “Users\Public\Public Documents\Thermo Fisher Scientific\AttuneNxt\Userdata” folder and exports the database to the specified location defined by the user.
- The database backup file that is created is called **attunenxt.backup** and includes information regarding the data structure. The data that are copied during the backup include copies of the Log files, Smart Monitor Data, and Userdata folders, which are located at the backup location.



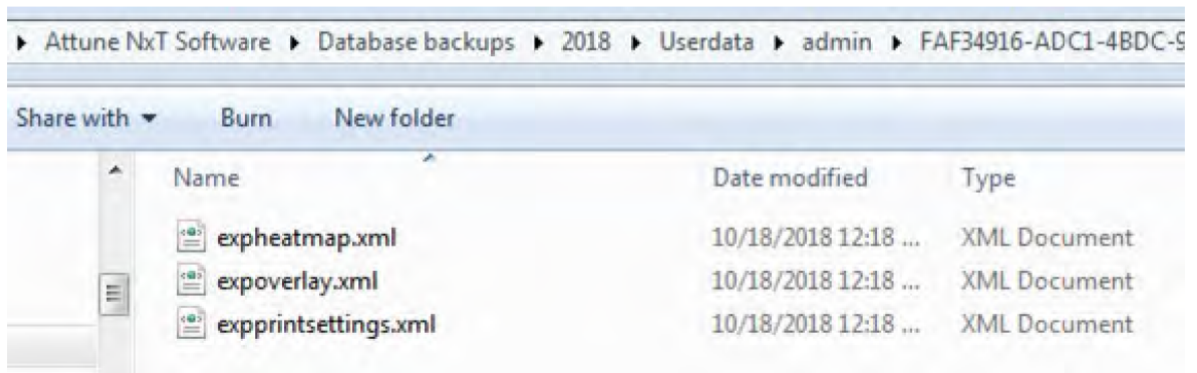
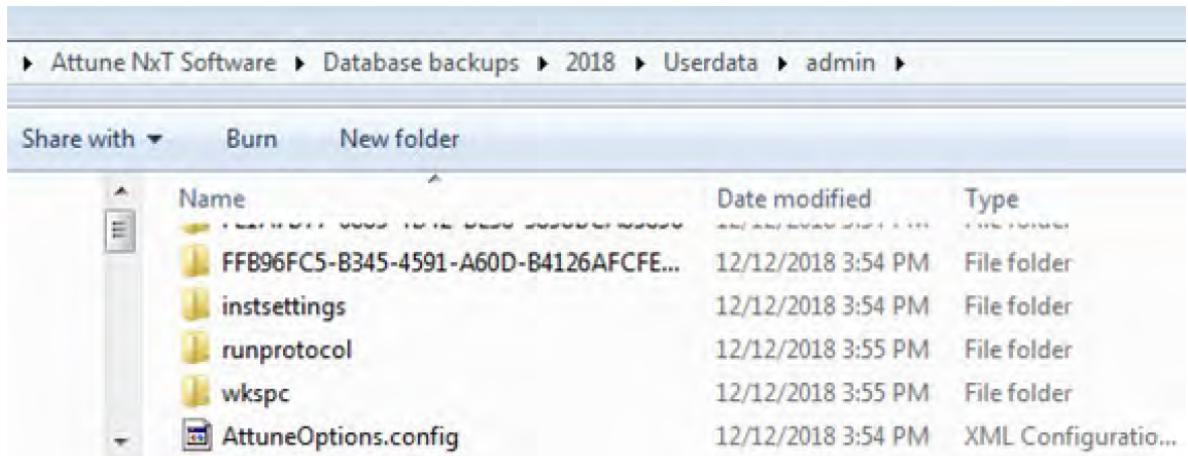
- Userdata folder contains subfolders for each account created in the software, including the default Service, System Administrator (sysadmin) and Administrator (admin) accounts, as well as custom, user-created accounts.





- In the folder for each user profile are data for the user account, including saved FCS files, and data associated with Experiment and Plate files, including Instrument settings, Workspaces, Compensation settings, Run protocols, Heat Map settings, Print settings, Overlay settings, Sample List settings, and Results Table settings.

The data are grouped into folders with easily identifiable names (Instrument settings, Run protocol settings, and Workspace settings) or saved in a folder with a Globally Unique Identifier label (GUID) that corresponds to the ID of the Sample or Experiment.





## Back up data

### Save and export Plate and Tube Experiment files

1. In the **Experiment Explorer**, right-click the **Experiment** or **Plate** name, then select **Export** ▶ **Experiment** (or **Plate**).

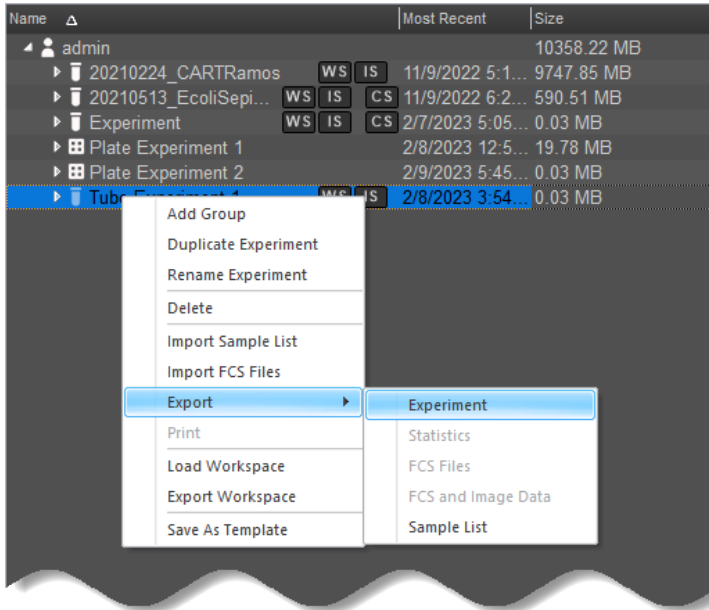


Figure 145 Export Tube Experiment from Experiment Explorer

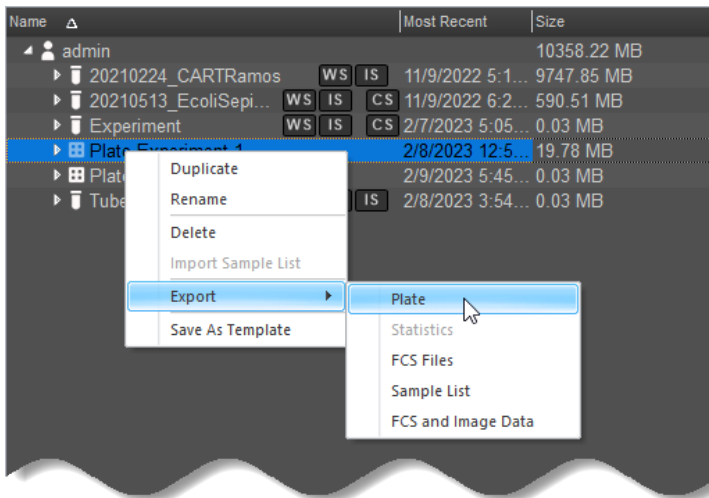
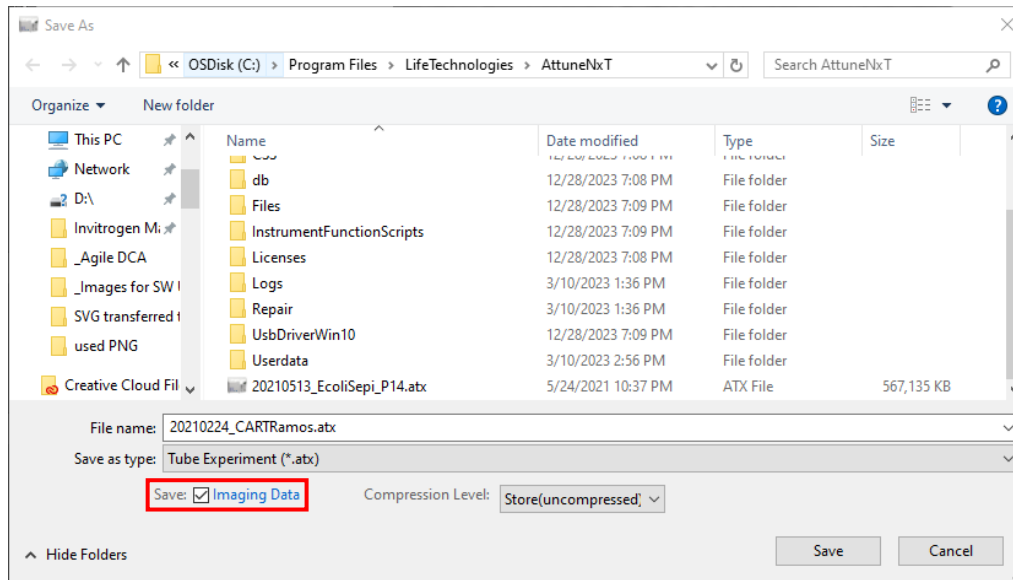


Figure 146 Export Plate Experiment from Experiment Explorer



2. In the **File Save (Export)** dialog, select the location to save the file.



3. To include image processing with the Experiment data that is being exported, ensure that the **Save Imaging Data** option is selected. To exclude the images and mask data from export, deselect the **Save Imaging Data** option.

---

**Note:** By default, the **Save Imaging Data** option is checked.

---

4. To save disk storage space, select **Compressed** from the **Compression Level** dropdown list.



---

**Note:** By default, **Store (uncompressed)** is selected for **Compression Level**.

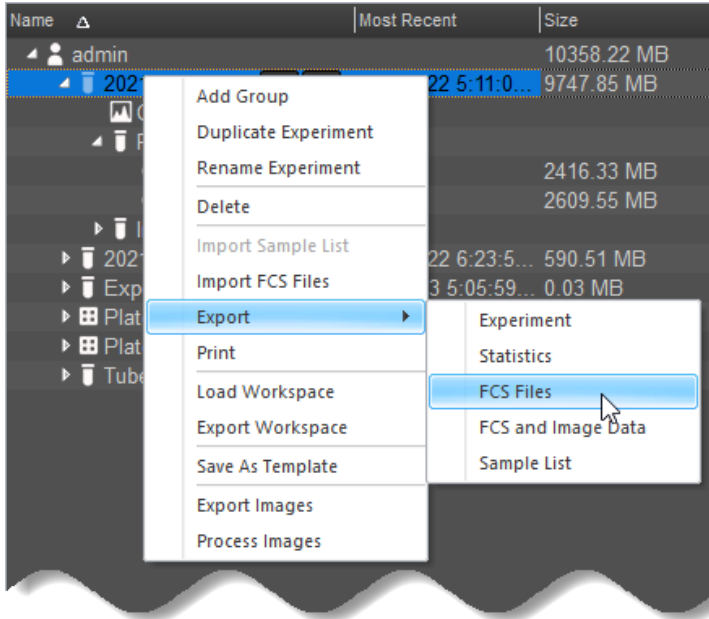
---

5. Click **Save** to save the file in the selected folder and close the dialog.

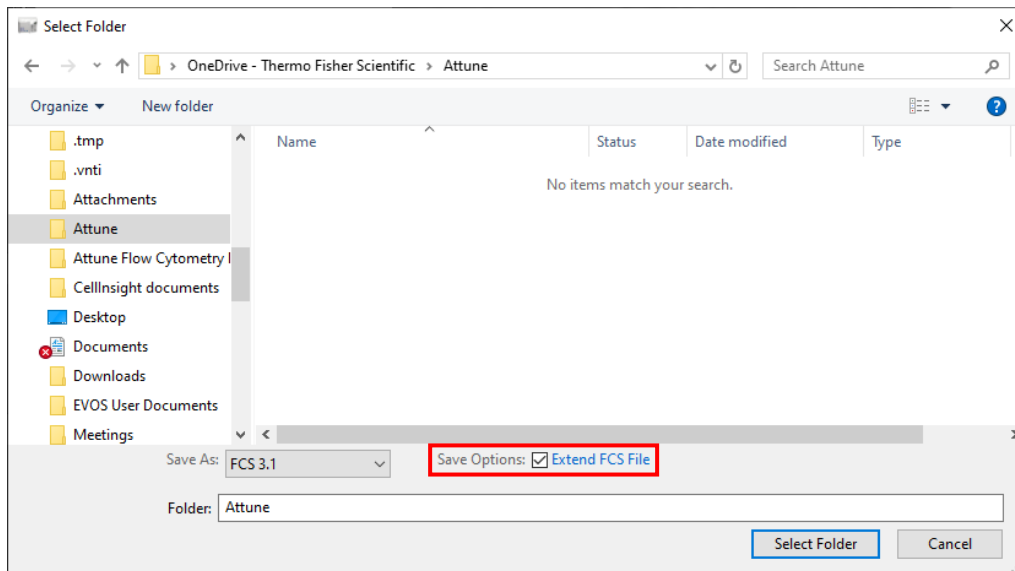


## Save and export FCS files

1. In the **Experiment Explorer**, right-click the **Experiment** or **Plate name**, then select **Export** ▶ **FCS Files**.



2. In the **Select Folder (Folder browser)** dialog, select the location to save the file, then click **Select Folder**.





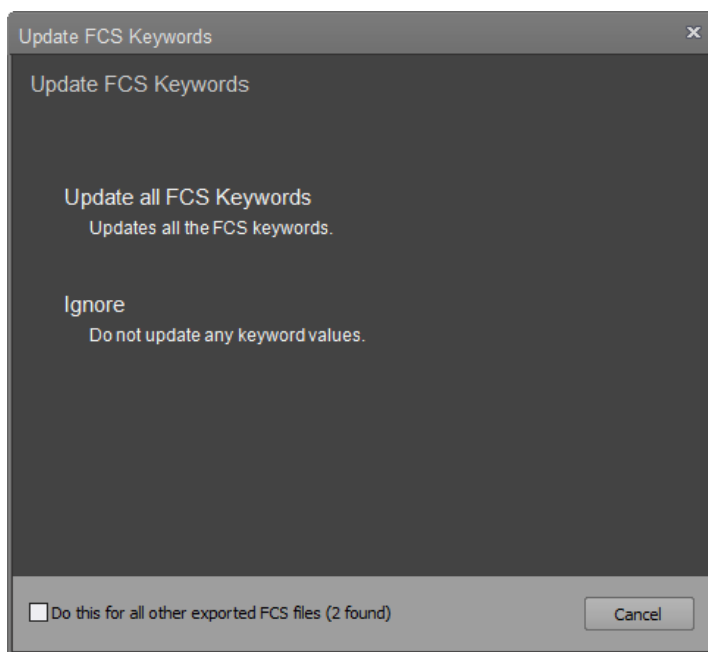
3. If desired, select **Extend FCS File** to add image processing results to the original FCS file, which creates a single FCS file that also includes the image processing parameters.

---

**Note:** When this option is selected, the original FCS file is appended with the image processing parameter data and keywords are updated so that the \$ORIGINALITY, \$LAST\_MODIFIER and \$LAST\_MODIFIED are set to indicate that the data have been appended and modified. The **Extend FCS File** option is available only if the Experiment has image processing data.

---

4. If desired, use the **Save As** dropdown change the FCS file type to **FCS 3.0**. By default, **FCS 3.1** is selected as the FCS file type.
5. In the **Update FCS Keywords** dialog, select **Update all FCS Keywords** to export FCS files with updated FCS keywords or select **Ignore** to export FCS files without updating any keyword values.



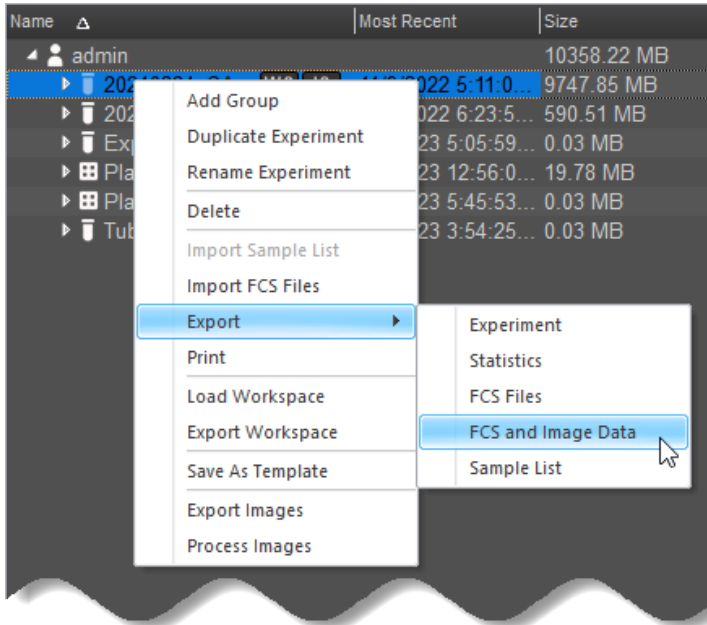
6. The software saves the FCS files in the selected folder, then closes the dialog.



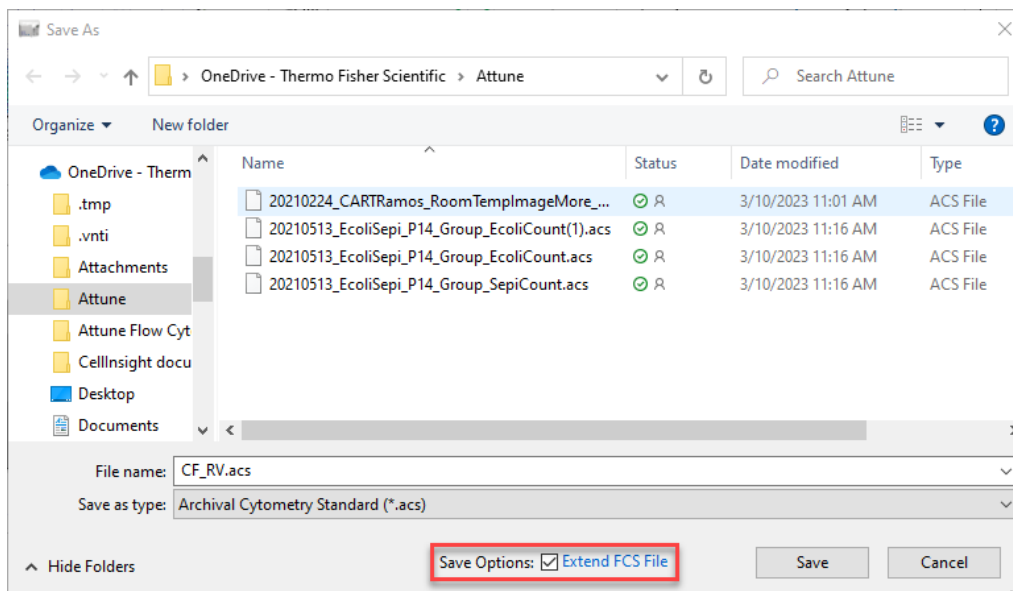
## Save and export FCS and image data

1. In the **Experiment Explorer**, right-click the **Experiment** or **Plate name**, then select **Export** ▶ **FCS and Image Data**.

**IMPORTANT!** The **Export** ▶ **FCS and Image Data** option is available only for experiments performed with an Attune™ CytPix™ Flow Cytometer, and it is active only if the experiment has FCS and image data.



2. In the **Folder Browser** dialog, select the location to save the file, then click **Select Folder**.







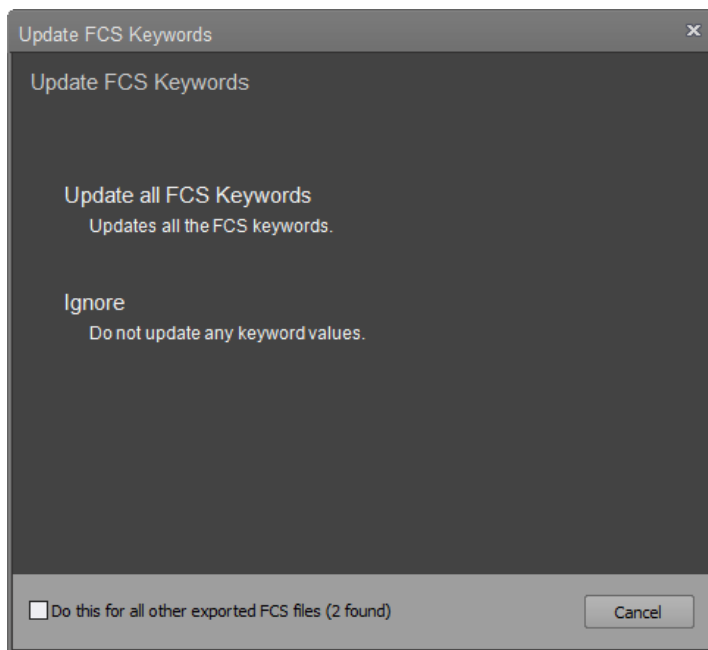
3. If desired, select **Extend FCS File** to add image processing results to the original FCS file, which creates a single FCS file that also includes the image processing parameters. This option is available only if the Experiment has image processing data.

---

**Note:** When this options is selected, the original the original FCS file is appended with the image processing parameter data and keywords are updated so that the \$ORIGINALITY, \$LAST\_MODIFIER and \$LAST\_MODIFIED are set to indicate that the data have been appended and modified.

---

4. In the **Update FCS Keywords** dialog, select **Update all FCS Keywords** to export FCS files with updated FCS keywords or select **Ignore** to export FCS files without updating any keyword values.



5. The software saves the FCS file, images, extended parameters (image processing data), and image mask data along with the FCS file into a single zipped file (based on ACS file format), then closes the dialog.

---

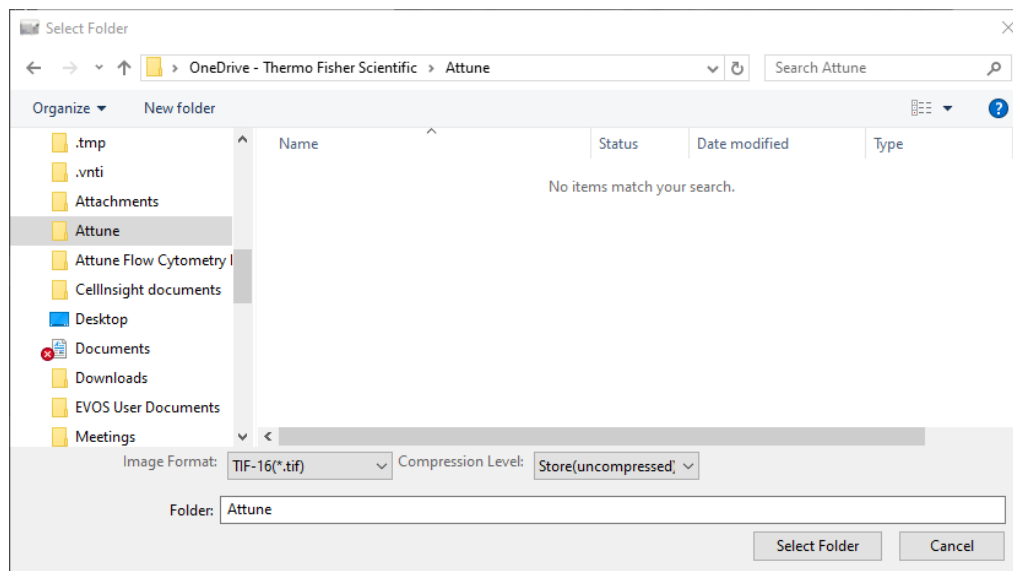
**Note:** Exported ACS files can be imported into a Sample, which imports the FCS data and all imaging data. Exported ACS files can be opened in third party software that supports the ACS standard (such as FlowJo™ Software and FCS Express™ Software).

---





2. In the **Select Folder (Folder browser)** dialog, select the location to save the file, then click **Select Folder**.



3. If desired, change the **Image Format** for the images to be exported.

---

**Note:** By default, images are exported as **TIF-16 (\*.tif)** files, which is the native file format collected from the Attune™ CytPix™ Flow Cytometer. You can select to save the images as **TIF-16, TIF-8, PNG, BMP, GIF, EMF, or JPG** files.

---

4. To save disk storage space, select **Compressed** from the **Compression Level** dropdown list.

---

**Note:** By default, **Store (uncompressed)** is selected for **Compression Level**.

---

5. The software saves the images from each Sample in the Experiment into a separate ZIP file in the selected folder, then closes the dialog.

---

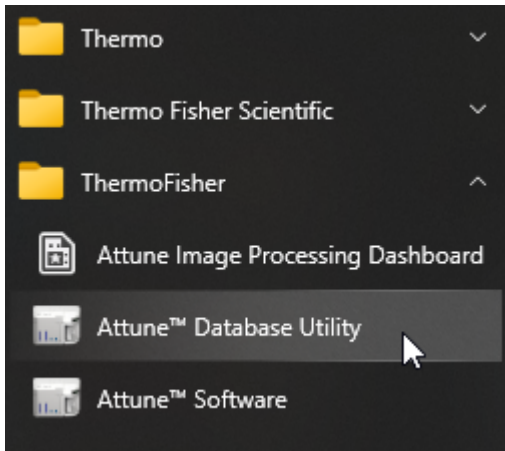
**IMPORTANT!** If images have been processed and mask overlays are turned on, images saved from the gallery or image view will have the mask data overlayed on the images.

---

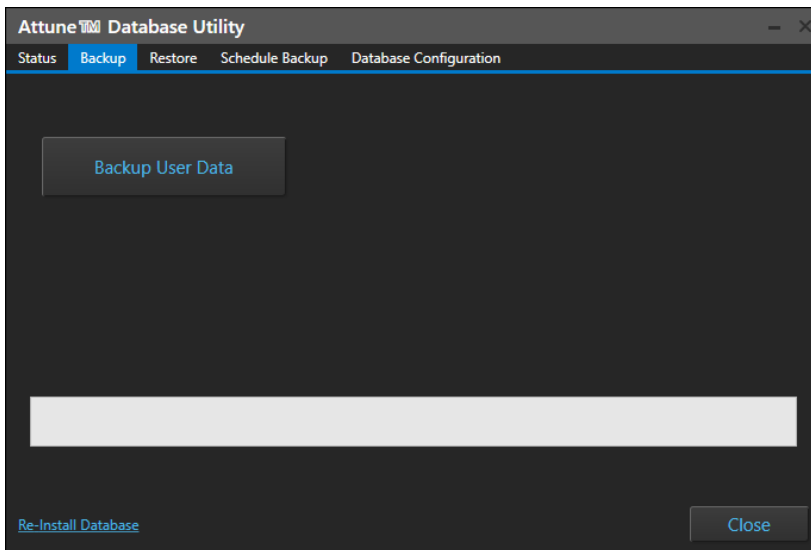


## Back up data using the Attune™ Database Utility

1. Close the Attune™ Cytometric Software.
2. Open the **Attune™ Database Utility** from the **Start Menu**, then sign in as **Administrator**.

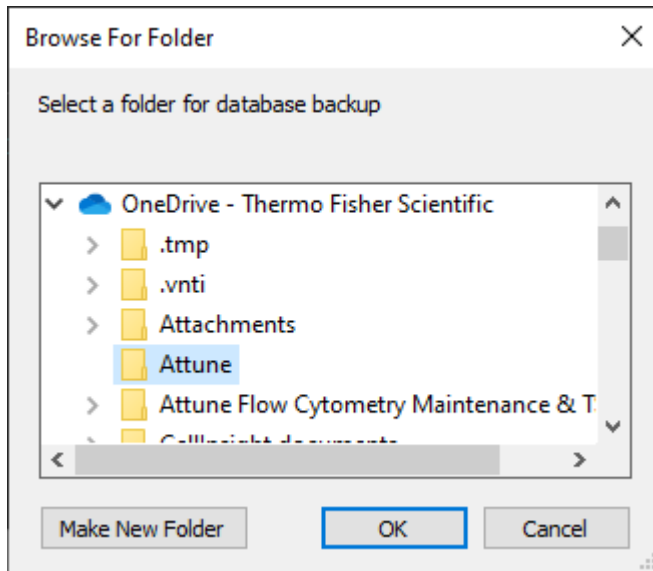


3. On the **Backup** tab, click **Backup User Data**.





4. Select the desired location to save the database.



As the backup continues, the progress bar indicates that files are being copied. When the backup is complete, the dialog states **Backup Complete**.



5. Click **Close** to exit the utility. A complete copy of all data associated with the Attune™ system has now been created.



## Restore data

### Restore data using the Attune™ Database Utility

You can use the Attune™ Database Utility to restore the software from a saved database. When you restore user data using a saved database, all data that are currently saved in the Attune™ Cytometric Software is replaced with data that were saved at the time the database backup was created. The database does not need to be created using the same version of software currently in use.

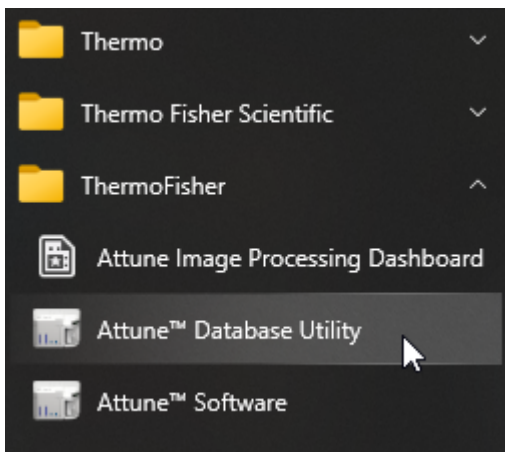
---

**IMPORTANT!** Restoration of user data from a saved database overwrites all data in the active, current instance of the Attune™ Cytometric Software.

Do **not** restore from a saved database without archiving data in the active instance of the Attune™ Cytometric Software.

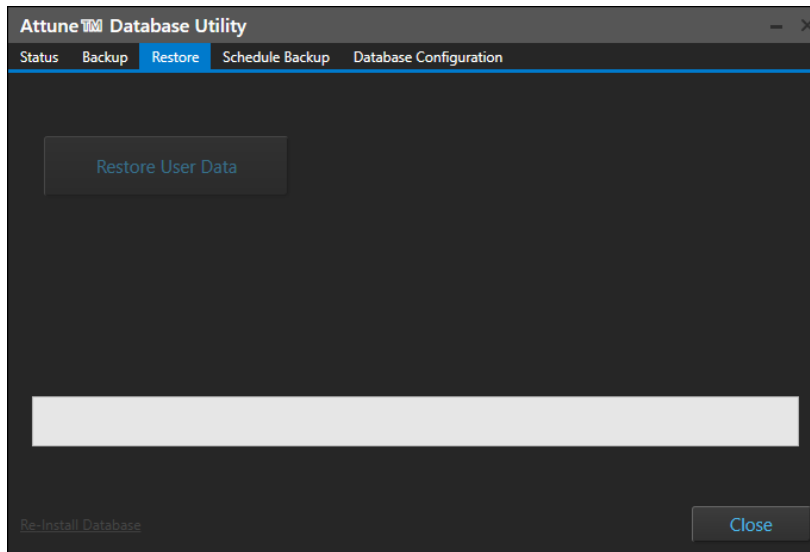
---

1. Close the Attune™ Cytometric Software.
2. Open the **Attune™ Database Utility** from the **Start** menu, then sign in as an **Administrator**.

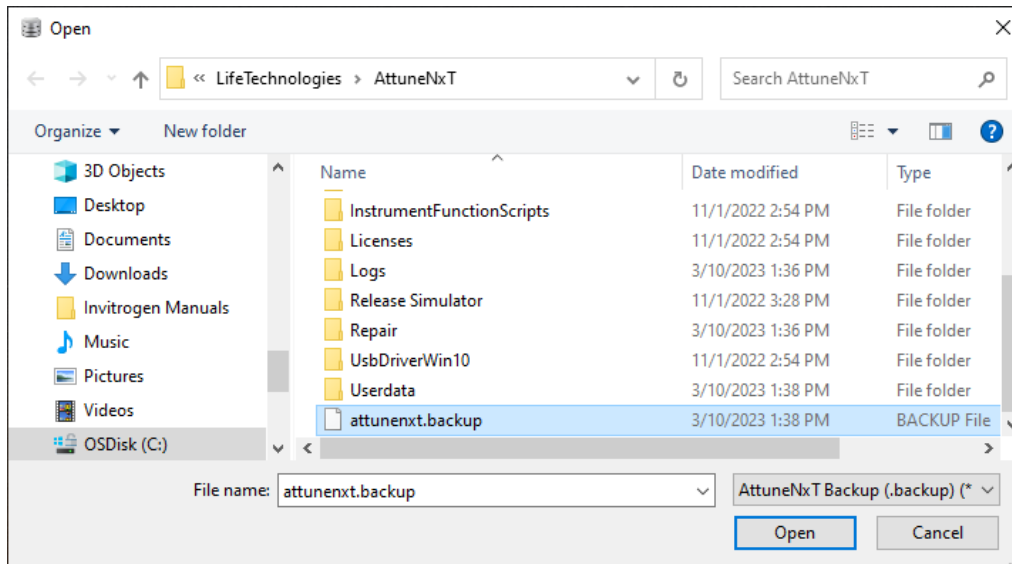




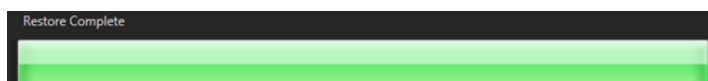
3. On the **Restore** tab, click **Restore User Data**.



4. In the **Open** dialog, select the correct database backup file (**attunenxt.backup**).



The database restoration begins immediately. As the restoration continues, the progress bar indicates the progress to completion. When the restoration is complete, the dialogue states **Restore Complete**.



5. Click **Close** to exit the utility. The Attune™ Cytometric Software now contains data, including Experiments, FCS files, user profiles, etc., from the database used for the restoration.



## Copy a database to a different computer

### Workflow

Copying an instance of the Attune™ Cytometric Software to a new computer is a 4-phase process that includes backing up data using the Attune™ Database Utility, copying all files to a new computer using an external hard drive or a network location, installing the Attune™ Cytometric Software on the new computer, and reinstalling the saved database.

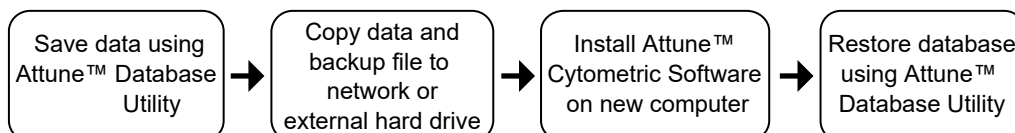
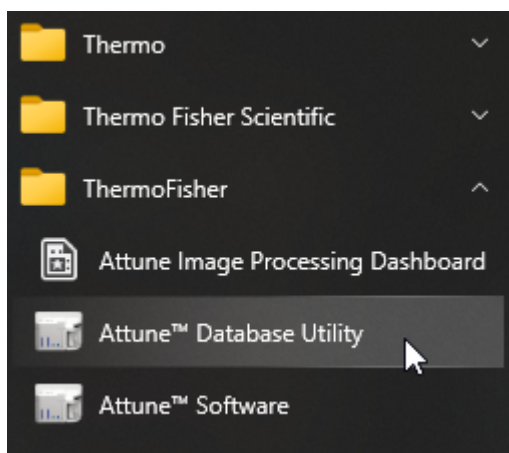


Figure 147 Workflow to copy an instance of the Attune™ Cytometric Software to another computer

### Save data using the Attune™ Database Utility

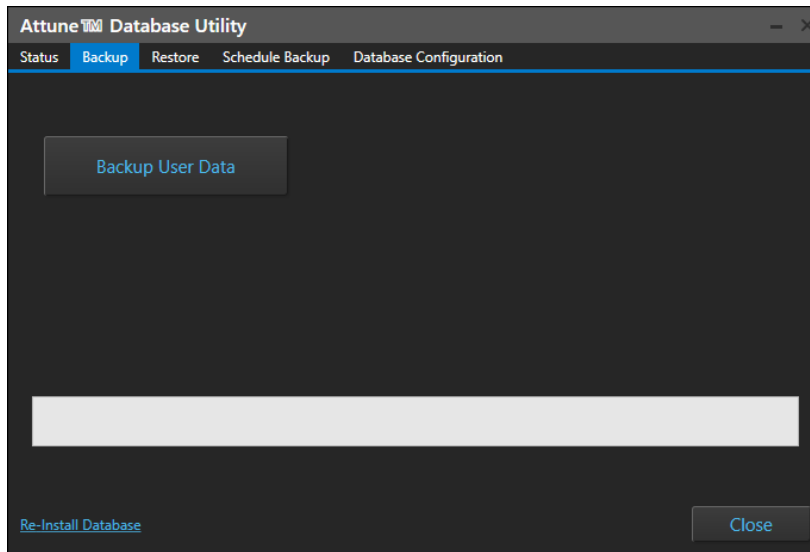
1. Close the Attune™ Cytometric Software.
2. Open the **Attune™ Database Utility** from the **Start** menu, then sign in as **Administrator**.



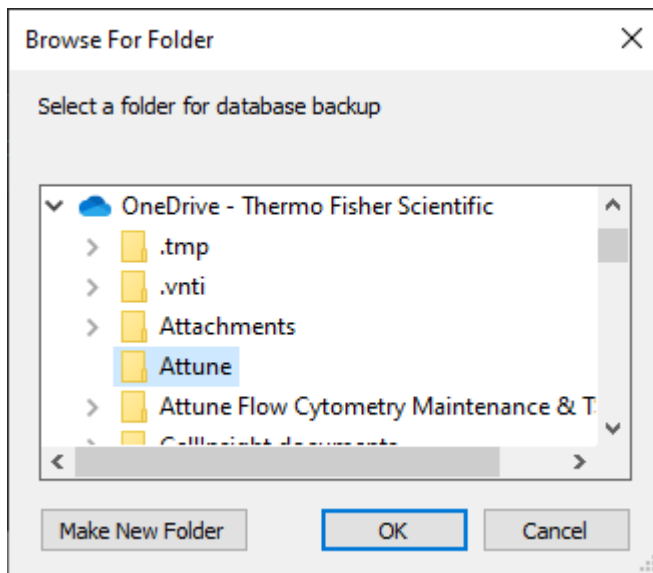




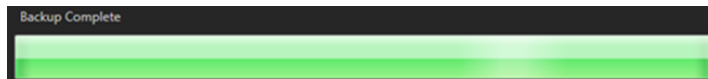
3. On the **Backup** tab, click **Backup User Data**.



4. Select the desired location to save the database.



As the backup continues, the progress bar indicates that files are being copied. When the backup is complete, the dialog states **Backup Complete**.

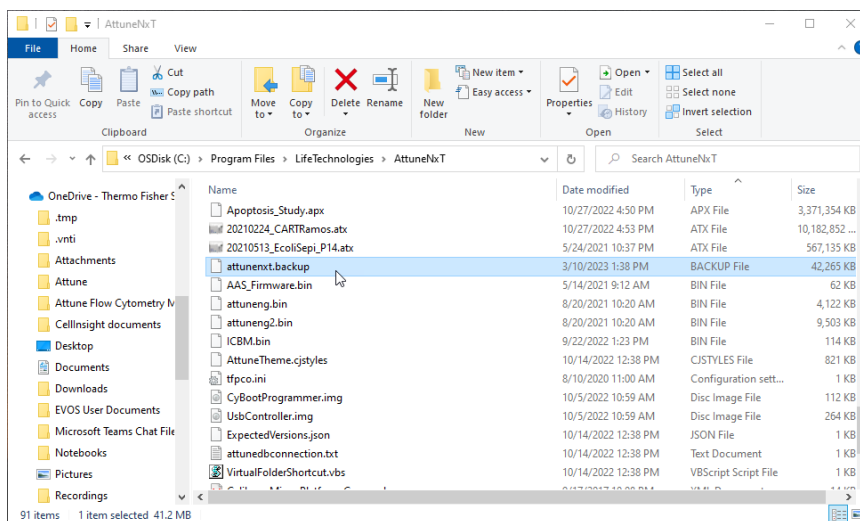


5. Click **Close** to exit the utility. A complete copy of all data associated with the Attune™ system has now been created.



## Copy data and backup file to network or external hard drive

1. Go to the location on the computer where the database file and associated data were saved.



2. Copy **all** data (folders and **attunenxt.backup** file) to another location (network location or external hard drive).

---

**IMPORTANT!** Failure to copy **all** data folders will cause database restoration to fail.

---

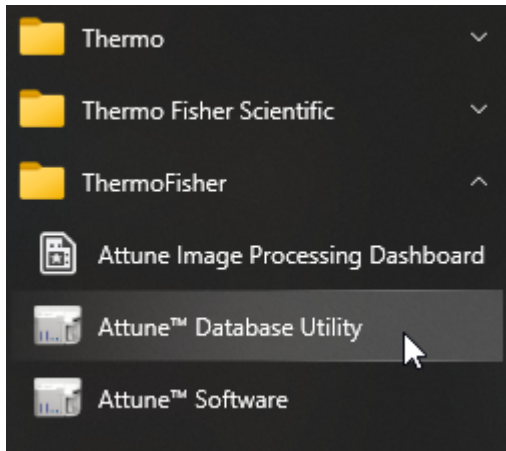
## Install Attune™ Cytometric Software on new computer

1. Install the Attune™ Cytometric Software on the secondary computer. For complete instructions, refer to the software release notes at [thermofisher.com/attune](https://thermofisher.com/attune).
2. If open, close the Attune™ Cytometric Software.

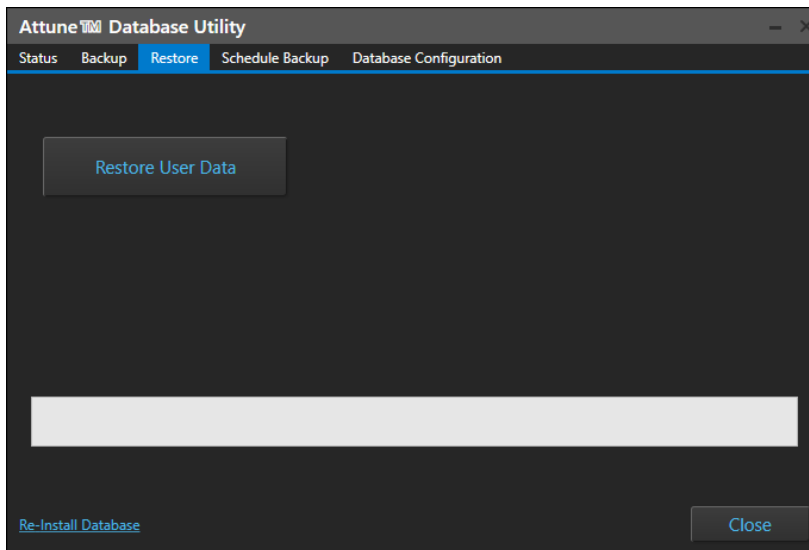


## Restore database using the Attune™ Database Utility

1. Copy all data folders and the **attunenxt.backup** file to the new computer.
2. Open the **Attune™ Database Utility** from **Start** menu, then sign in as an **Administrator**.

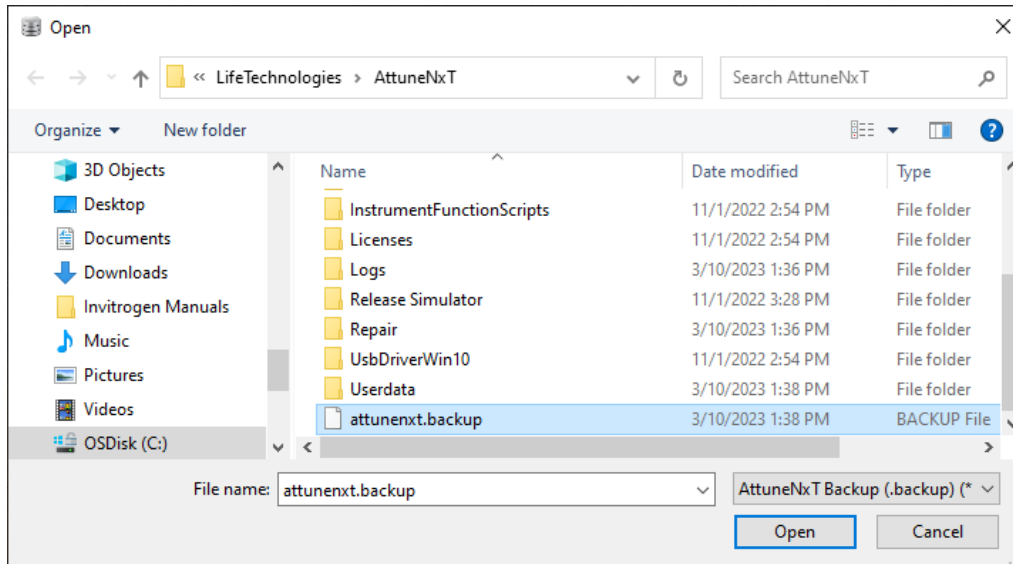


3. On the **Restore** tab, click **Restore User Data**.





4. In the **Open** dialog, select the database backup file (**attunenxt.backup**) that was saved on the computer.



The database restoration begins immediately. As the restoration continues, the progress bar indicates the progress to completion. When the restoration is complete, the dialogue states **Restore Complete**.

5. Click **Close** to exit the utility. The Attune™ Cytometric Software Cytometric Software now contains data, including Experiments, FCS files, user profiles, etc., from the database used for restoration.



# Permanently remove saved data from the Attune™ computer

## Data deletion options

The recommended method for deletion of data from the Attune™ Cytometric Software is to delete Tube or Plate Experiments from the Experiment Explorer in each user profile. When deleting files from individual user accounts is not possible, you can use one the other options described in this section.

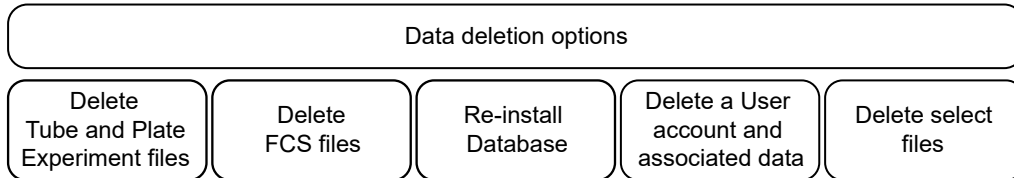


Figure 148 Options for permanently removing saved data from the Attune™ computer

## Delete Tube and Plate Experiment files

1. Sign in to the account that includes the data you want to delete.
2. Right-click the **Experiment**, **Plate**, or **Sample** name, then select **Delete** from the context menu.

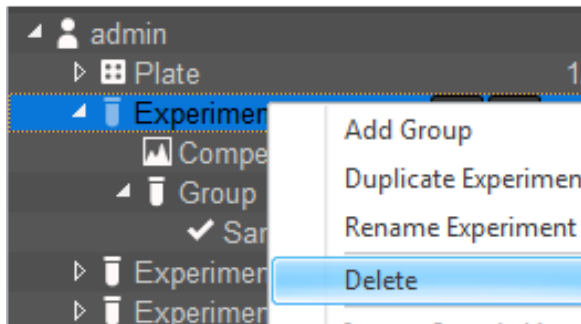


Figure 149 Delete Tube Experiment

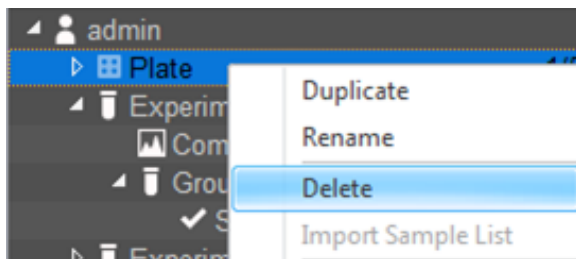


Figure 150 Delete Plate Experiment

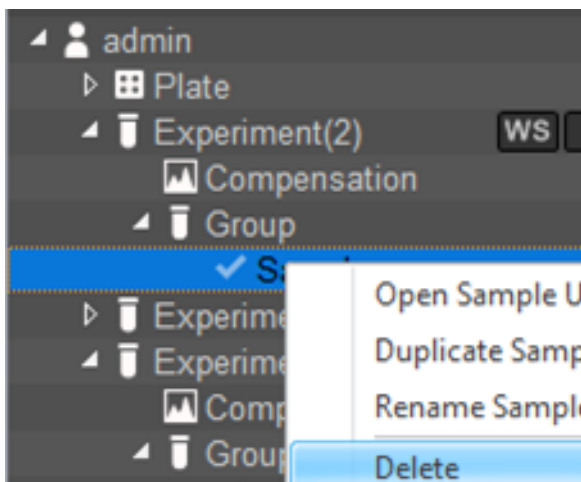


Figure 151 Delete Sample

3. *Optional:* To select multiple Tube Experiments, Plate Experiments, or FCS files for bulk deletion, select each file individually as you hold the keyboard **Ctrl** key.

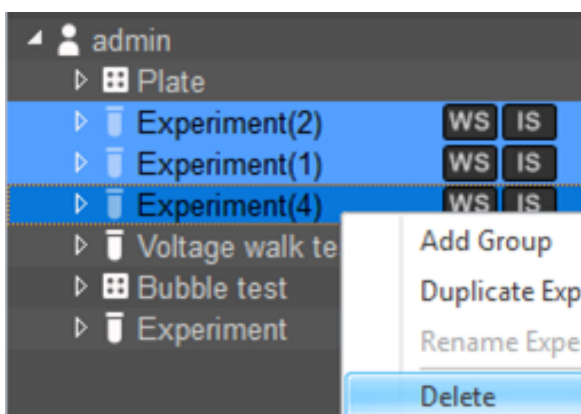


Figure 152 Delete multiple Tube Experiments

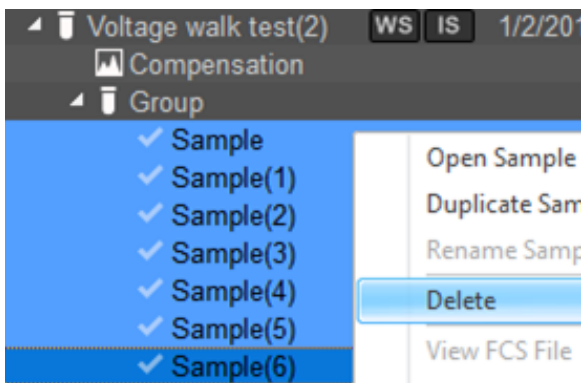
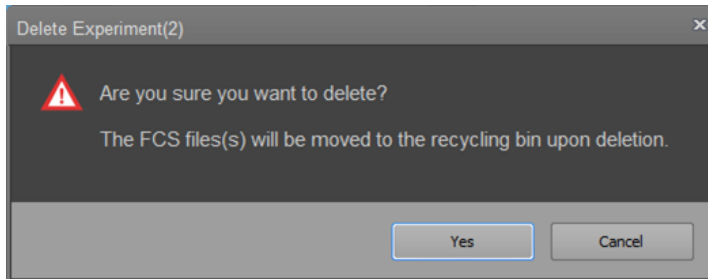


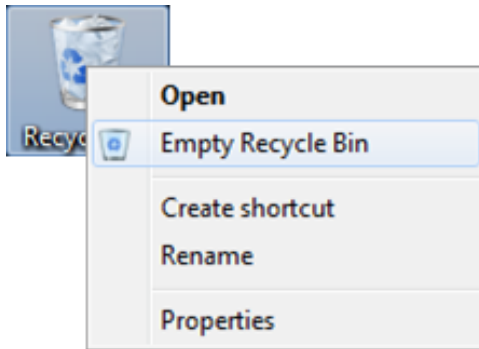
Figure 153 Delete multiple Samples



4. In the Delete dialog, click **Yes** to confirm the file deletion or click **Cancel** to close the dialog without deleting the selected files. The files that are deleted from the software are sent to the **Recycle Bin** of the computer.



5. Empty the **Recycle Bin** to remove the deleted files permanently from the computer. Alternatively, open the **Recycle Bin** and select individual files to delete permanently.





## Reinstall a new database to erase existing data

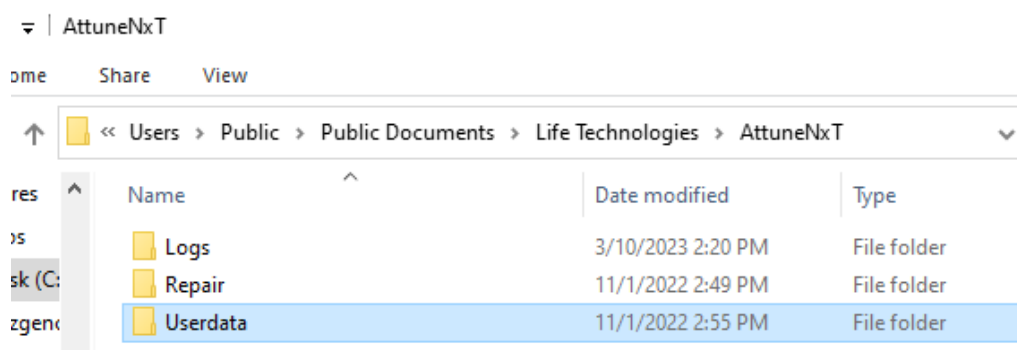
You can use the Attune™ Database Utility to reinstall a new, empty database. Installation of a new database completely erases all existing data and user accounts in the current, active instance of the Attune™ Cytometric Software.

---

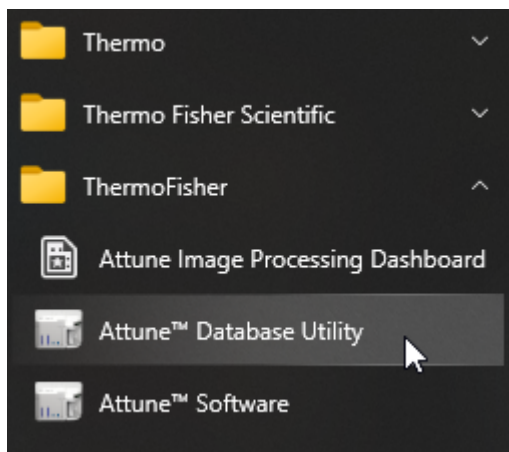
**IMPORTANT!** Do **not** reinstall a new database without first saving and exporting the experiments and files from the Attune™ Cytometric Software.

---

1. Close the Attune™ Cytometric Software.
2. Go to the **C:\Users\Public\Public Documents\Life Technologies\AttuneNxT** folder, then right-click the **Userdata** folder and select **Delete** from the menu.



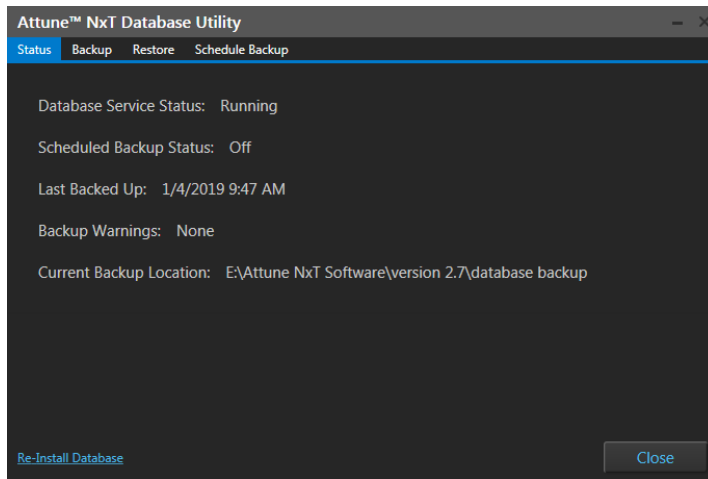
3. Open the **Attune™ Database Utility**, then sign in as an **Administrator**.



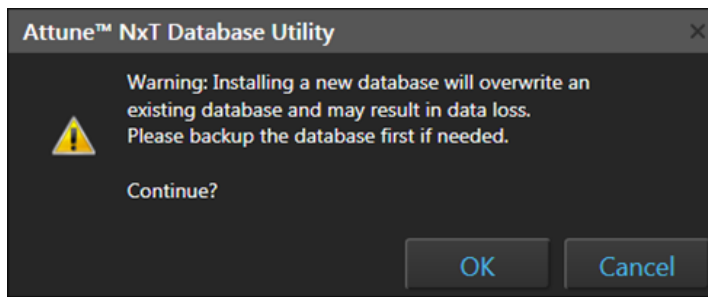




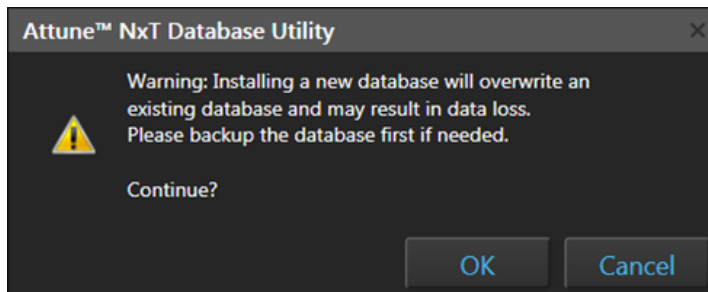
4. In the **Status** tab, select **Re-install Database** (in the lower left corner).



5. Click **OK** to confirm the intent to install a new database or click **Cancel** to close the dialog without installing a new database.



6. When the new database installation is completed, click **Close**.



7. Open the Attune™ Cytometric Software. The new database does not include any data from the previously created accounts.

---

**Note:** Sign in to the Attune™ Cytometric Software using one of the default account settings (**sysadmin** or **admin**).

The default username for the Administrator account is **admin** and the default password is **admin**.

The default username for the System Administrator account is **sysadmin** and the default password is **sysadmin**.

---



## Delete a user account to remove data

The recommended method for deletion of data from the Attune™ Cytometric Software is to delete Tube or Plate Experiments from the Experiment Explorer in each user profile. However, if a user account is no longer needed, you can delete the account to remove all files associated with it.

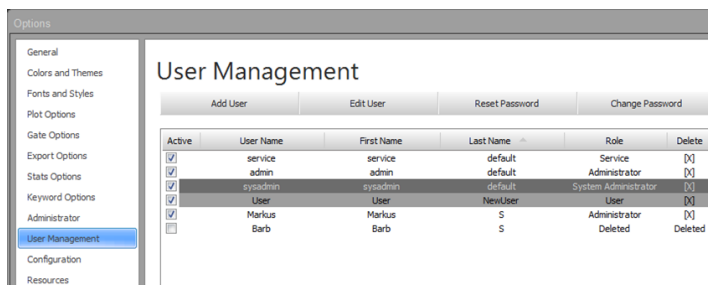
**IMPORTANT!** Deleting a user account permanently removes the user account and all associated files from the Attune™ Cytometric Software.

We do **not** recommend deletion of a user account without first backing up the user data to an external location.

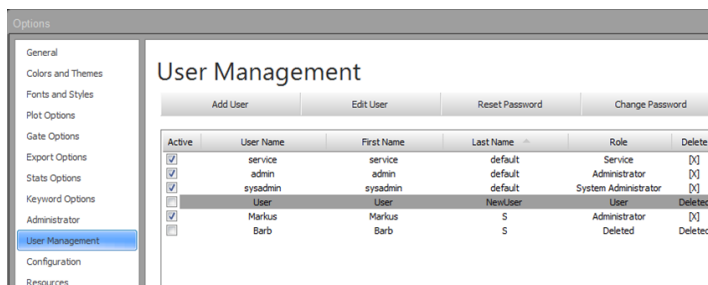
1. Sign in to the Attune™ Cytometric Software as the **System Administrator (sysadmin)**.

**Note:** The default username for the System Administrator account is **sysadmin** and the default password is **sysadmin**.

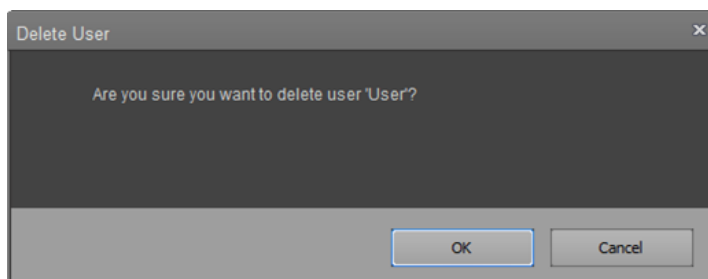
2. In the **Options** dialog, select **User Management**.



3. Select the user account you want to delete, then click the **X** in the **Delete** column.

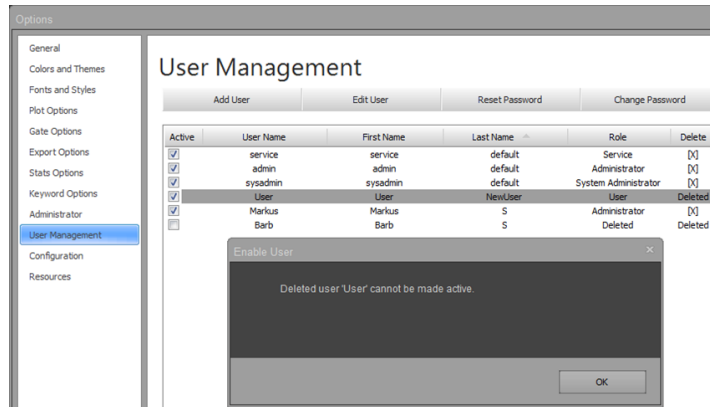


4. In the **Delete User** dialog, click **OK** to confirm the intent to delete the selected user account or click **Cancel** to close the dialog without deleting the account.





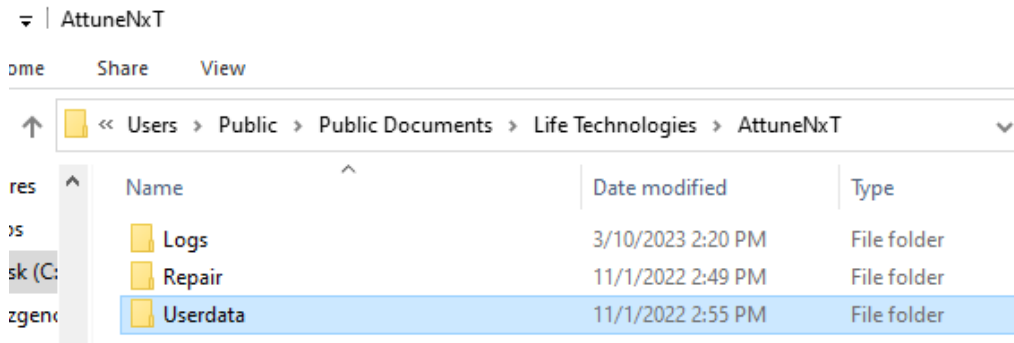
When the account has been deleted, the Active box will be unchecked and the user status cannot be reenabled by selecting the Active box.



- Although the user account has been deleted, all associated user data remains in the Attune™ NxT database until the data are manually deleted.

To delete all associated data, go to

**C:\Users\Public\Documents\Life Technologies\AttuneNxT\Userdata\X**, where X is the user account name for the deleted user account.



- Right-click the data folder for the deleted user account, then select **Delete** from the dropdown menu.
- In the **Delete Folder** dialog, click **Yes** to confirm the intent to move the data folder to the **Recycle Bin** or click **No** to close the dialog without deleting.
- Empty the **Recycle Bin** to remove the deleted files permanently from the computer. Alternatively, open the **Recycle Bin** and select individual files to delete permanently.

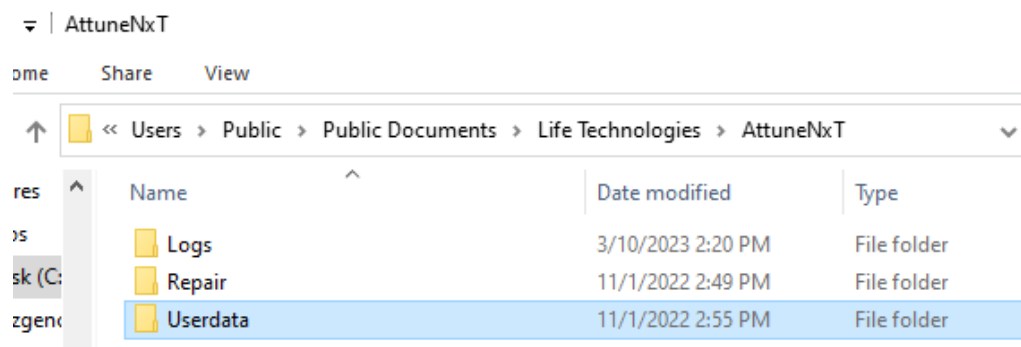
---

**Note:** When a user account is deleted, the folder corresponding to the deleted account continues to be visible in C:\Users\Public\Documents\Life Technologies\AttuneNxT\Userdata. This folder can be opened, but it contains empty subfolders and no actual data.



## Appendix C Data management in Attune™ Cytometric Software

Permanently remove saved data from the Attune™ computer





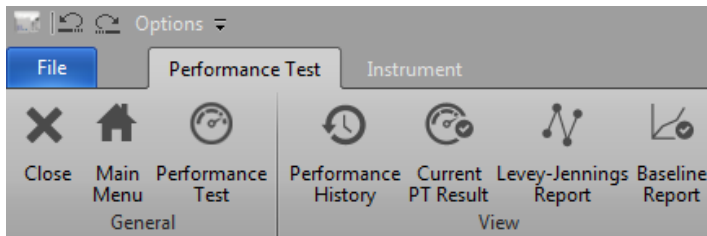
# Permanently remove old baseline and performance test data

## About baseline and performance test data

Baseline and Performance Test generate many data files including XML files associated with laser delay and test results, FCS files, CSV files, and log files annotating steps that occur during Baseline and Performance Test.

BL-2019-02-13T16.19.18.000	2/13/2019 4:22 PM	File folder
BL-2019-02-13T16.33.44.000	2/13/2019 4:37 PM	File folder
PT-2019-02-13T15.53.01.000	2/13/2019 3:54 PM	File folder
PT-2019-02-13T15.55.00.000	2/13/2019 3:56 PM	File folder
PT-2019-02-13T16.03.13.000	2/13/2019 4:04 PM	File folder
BL-2019-02-13T14.55.43.000LaserDelay.xml	2/13/2019 3:51 PM	XML File
BL-2019-02-13T14.55.43.000Results.xml	2/13/2019 3:51 PM	XML File
BL-2019-02-13T16.19.18.000Results.xml	2/13/2019 4:22 PM	XML File

Baseline and Performance Test data are stored in two locations in the Attune™ Cytometric Software: directly in the database and in the Public Documents folder. In the Attune™ Cytometric Software, data are pulled directly from the database to view Baseline or Performance Test Results, Levey-Jennings data, Performance History, and the Baseline Report.





# Appendix C Data management in Attune™ Cytometric Software

Permanently remove old baseline and performance test data

Options <admin> - Attune™ NxT Software v3.1.0

File Performance Test Instrument

Close Main Menu Performance Test Performance History Current PT Result Levey-Jennings Report Baseline Report

## Baseline Test Results

Baseline test successful

Baseline: 756080D - 2/13/2019 2/13/2019 2:55:43 PM

Channel	PMTV	Target MFI	MFI	Robust %CV	Qr	Background	Linearity	ASF	Laser Delay	Result
FSC	487	300,000	298,796	1.68 %	0.000	0	0.000	1.11	1100	✓
SSC	285	300,000	294,170	3.21 %	0.000	0	0.000	1.11	1100	✓
BL1	378	300,000	295,691	1.24 %	0.047	127	1.000	1.11	1100	✓
BL2	311	300,000	295,495	1.17 %	0.065	169	0.966	1.11	1100	✓
BL3	349	300,000	299,207	1.82 %	0.049	39	1.000	1.11	1100	✓
RL1	333	300,000	295,093	3.84 %	0.066	100	0.997	1.04	1497	✓
RL2	353	300,000	296,932	3.60 %	0.010	306	1.000	1.04	1497	✓
RL3	423	300,000	297,819	4.01 %	0.047	141	0.998	1.04	1497	✓
VL1	297	300,000	300,000	2.80 %	0.010	2738	1.000	1.14	702	✓
VL2	325	300,000	294,262	2.64 %	0.010	1001	0.997	1.14	702	✓
VL3	324	300,000	296,212	2.78 %	0.016	182	0.996	1.14	702	✓
VL4	397	300,000	297,477	3.42 %	0.003	298	0.987	1.14	702	✓
YL1	381	300,000	294,021	1.33 %	0.081	37	0.999	1.12	305	✓
YL2	348	300,000	297,493	1.27 %	0.046	68	0.972	1.12	305	✓
YL3	392	300,000	299,462	2.07 %	0.035	134	0.999	1.12	305	✓
YL4	400	300,000	296,013	2.58 %	0.006	160	1.000	1.12	305	✓

## Levey-Jennings Report

Baseline: 1759476A - 8/7/2017 8/7/2017 9/1/2017

Plotting Statistic: PMT voltage Laser Scatter Blue Red Violet Yellow

### FSC PMT voltage

Date	PMTV
7 Aug	380
10 Aug	380
13 Aug	380
16 Aug	380
19 Aug	380
22 Aug	380
25 Aug	380
28 Aug	380
31 Aug	380

August 2017 - September 2017

### SSC PMT voltage

Date	PMTV
7 Aug	320
10 Aug	320
13 Aug	320
16 Aug	320
19 Aug	320
22 Aug	320
25 Aug	320
28 Aug	320
31 Aug	320

August 2017 - September 2017

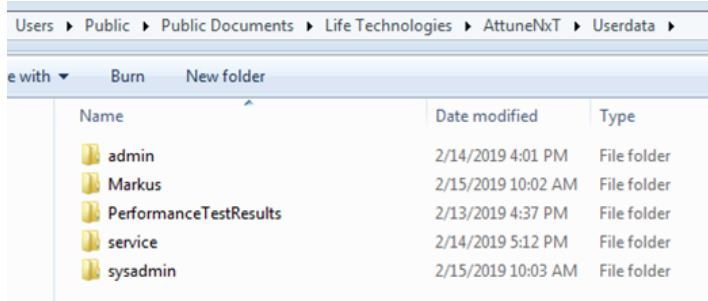


Over time, the PerformanceTestResults folder (X:\Users\Public\Documents\Life Technologies\AttuneNxT\Userdata) can become very large. To permanently remove the files contained in the PerformanceTestResults folder, follow the steps described in this section.

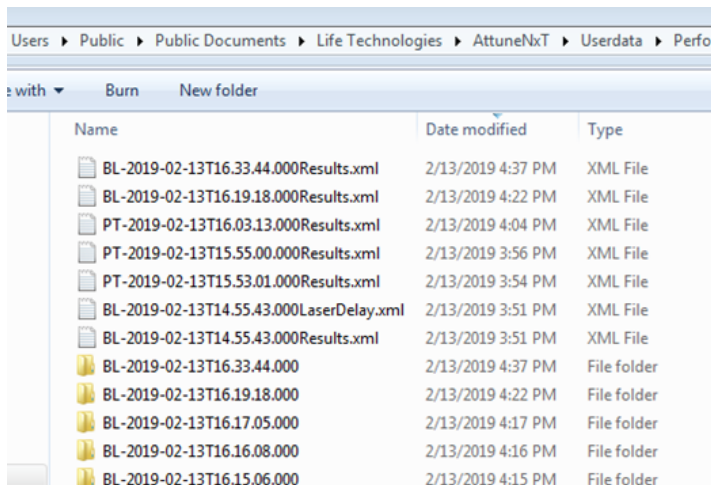


## Remove baseline and performance test data in Attune™ Cytometric Software version 2.7 or earlier

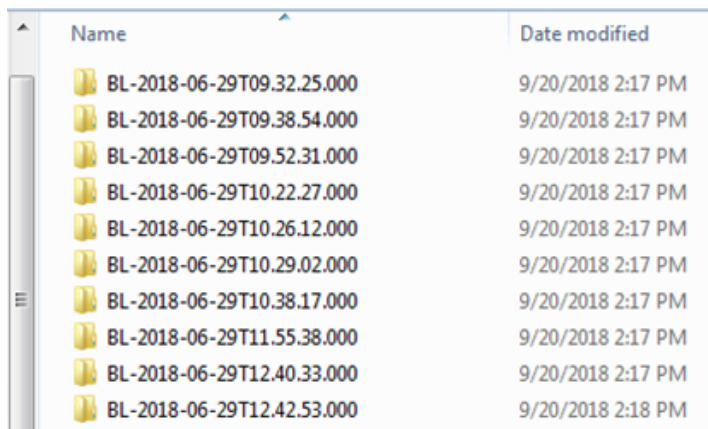
1. Go to the **PerformanceTestResults** folder  
(X:\Users\Public\Documents\Life Technologies\AttuneNxT\Userdata).



2. In the PerformanceTestResults folder, organize data based on the “Date modified column” by clicking on the **Date modified column header**.



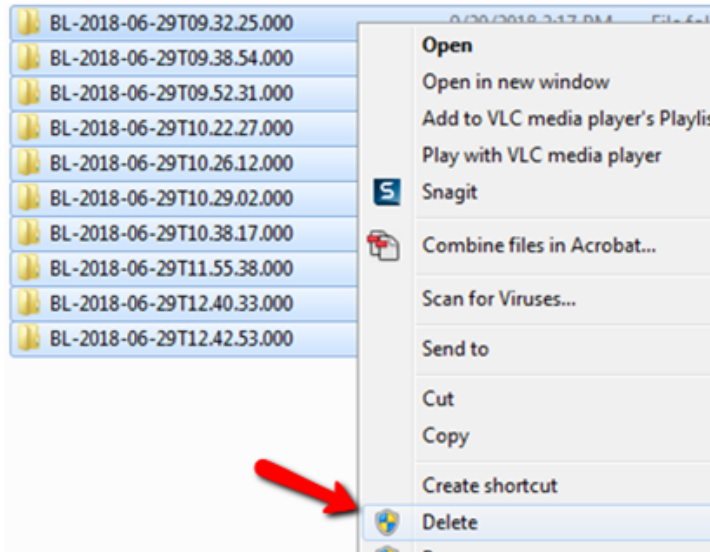
3. Select data to delete based on the date the data was created. We recommend keeping the last 3 months of data in the folder in case these files are needed for troubleshooting. In the following example, select the files created before 13 November 2018.



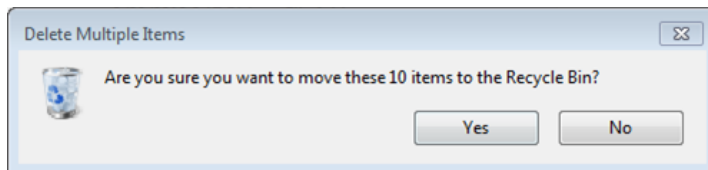




- Multiselect the appropriate files to delete, right-click, then select **Delete** from the dropdown menu to delete the files.



- Click **Yes** to confirm the deletion.

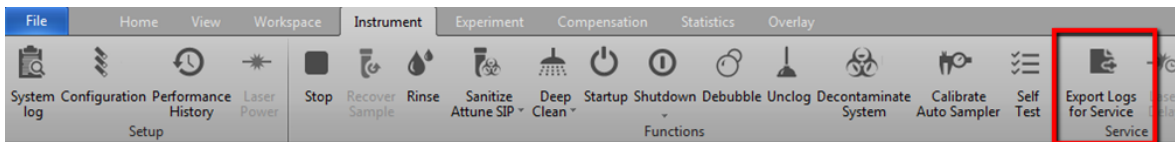


- Empty the **Recycle Bin** of the computer to remove the deleted files permanently from the computer.

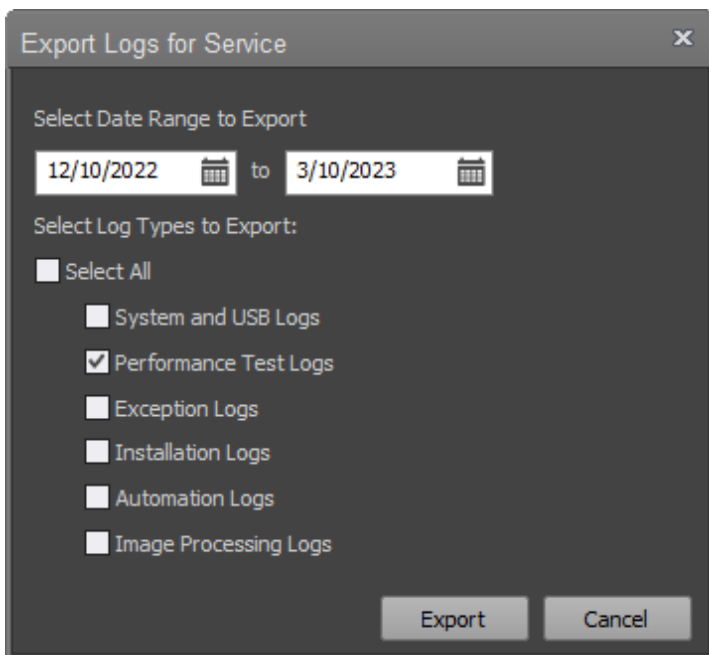


## Remove baseline and performance test data in Attune™ Cytometric Software version 3.1 and later

1. Open the Attune™ Cytometric Software and sign in to an administrator account.
2. Archive at least 3 months of Baseline and Performance Test Data before deletion. To do this, click **Export Logs for Service** on the **Instrument** tab of the **Ribbon bar**.

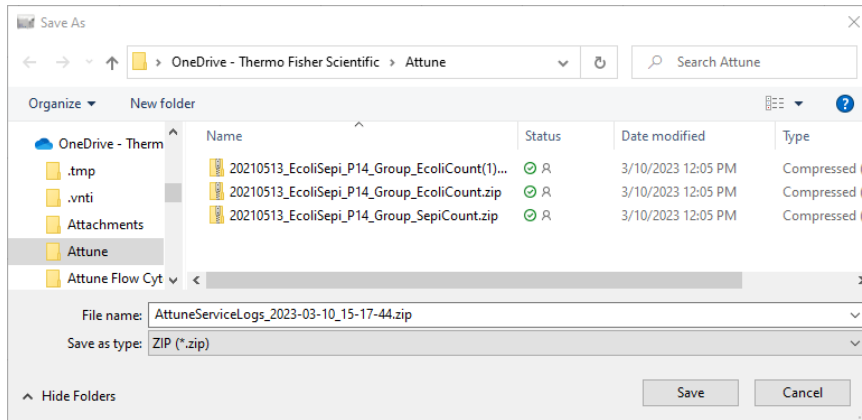


3. In the **Export Logs for Service** dialogue, select **Date Range to Export**, check the **Performance Test Logs** option, then click **Export**.

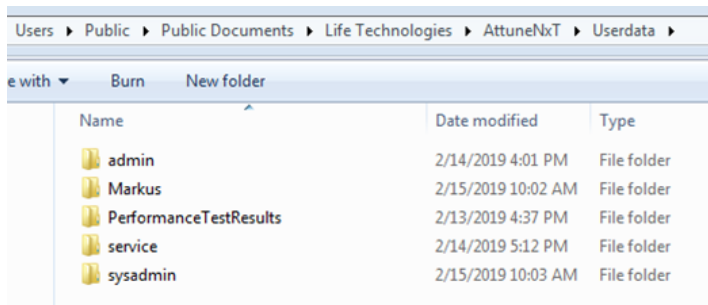




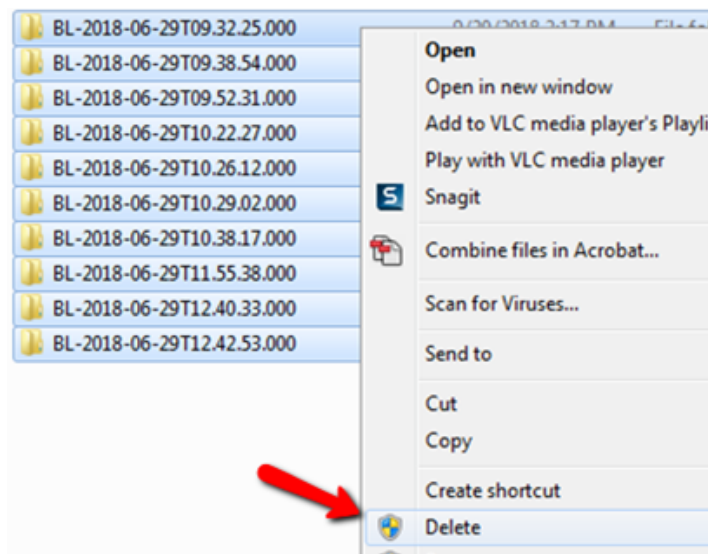
4. In the **Save As** dialogue, select the file to save the data, then click **Save**. All files required for troubleshooting are exported to the save location as a zip folder.



5. Go to the **PerformanceTestResults** folder (X:\Users\Public\Documents\Life Technologies\AttuneNxT\Userdata).

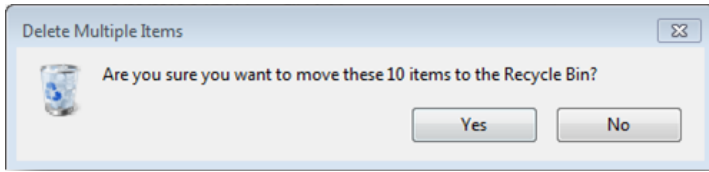


6. Open the **PerformanceTestResults** folder, multiselect all files in the folder, right-click, then select **Delete** from the dropdown menu to delete the files.





7. Click **Yes** to confirm the deletion.



8. Empty the **Recycle Bin** of the computer to remove the deleted files permanently from the computer.



# SAE Administrator Console

## SAE module

### Overview of the SAE module

The **SAE module** (Security, Audit, and Electronic Signature) is a client-server software configuration that is used to meet specific requirements for security, audit, and e-Signature (such as 21 CFR Part 11 compliance).

---

**IMPORTANT!** 21 CFR part 11 is a regulation that describes the criteria for acceptance by the FDA for electronic records and electronic signatures. Part 11 is composed of procedural and technical requirements. Procedural requirements are the standard operating procedures instituted by the end user, and technical requirements are the technical characteristics of the software used.

The **SAE module** of the Attune™ Cytometric Software does not automatically guarantee 21 CFR part 11 compliance. Compliance is the consequence of the **end user's work process and systems used**. Attune™ Cytometric Software in the **SAE mode** enables 21 CFR part 11 compliance for the flow cytometry data collection and analysis steps in the workflow.

---

The **SAE module** includes three components:

- **SAE Administrator Console**—Tool that is used by an SAE administrator to configure the **SAE module** (see “SAE Administrator Console” on page 880).
- **SAE server (server)**—Service that runs in the background and stores SAE settings, user accounts, audit records, and e-Signature records. By default, the SAE server is installed on the same computer as the SAE Administrator Console.
- **SAE screens (client)**—Screens that are displayed (sign in, audit, and e-Signature) in the application (such as the Attune™ Cytometric Software) and that require user input.

The **SAE module** can be configured to provide the following functionality:

Function	Description
System Security	Controls user access to the Attune™ application. A default user account with the <b>Administrator</b> role is provided at installation that has access to both SAE Administrator Console and the Attune™ Cytometric Software. Other default SAE user roles include <b>Advanced User, User, Reviewer, and No Privileges</b> roles (“SAE account permissions” on page 887). You can set up additional SAE user accounts with specific permissions.

(continued)

Function	Description
Auditing	Tracks actions performed by users, changes to SAE settings, and changes made to application objects (i.e., Experiments) (“Audit tab” on page 903). You can: <ul style="list-style-type: none"> <li>• Enable or disable audits, select actions and application objects to be audited, specify the audit mode, and require users to provide a reason for changes made to auditable application objects.</li> <li>• View specific audit logs and generate printable audit records.</li> <li>• Archive the audit records or configure auto archive settings.</li> </ul>
Electronic Signature (e-Signature)	Determines if users are required to fulfill signature requirements before performing specific functions. You can: <ul style="list-style-type: none"> <li>• Configure e-signature so that a user can export data, print data, and save experiment as a template only if the required e-Signatures are provided.</li> <li>• Configure each e-Signature event to require multiple signatures and to require users with specific roles to sign.</li> </ul>

## SAE Administrator Console

### Overview of the SAE Administrator Console

The **SAE Administrator Console** is used to configure the **SAE module**, including user credential management (username and password complexity), users, user roles (security), auditing, and e-Signature features.

In the SAE Administrator Console, a software or instrument that is configured for the **SAE module** is called an “**application**”. An example application is the Attune™ Cytometric Software.

The SAE Administrator Console has an Attune™-specific security profile that defines application permissions, default user roles, auditable application objects, signature actions, signature application objects, and default signature meanings.

Changes made on the SAE Administrator Console are synchronized with the Attune™ Cytometric Software every 10 seconds. If the SAE Administrator Console is offline, then the Attune™ Cytometric Software uses a cache to manage a local copy of the SAE Administrator Console server settings and a cache of all audit and e-Signature records that are created while the server is offline.

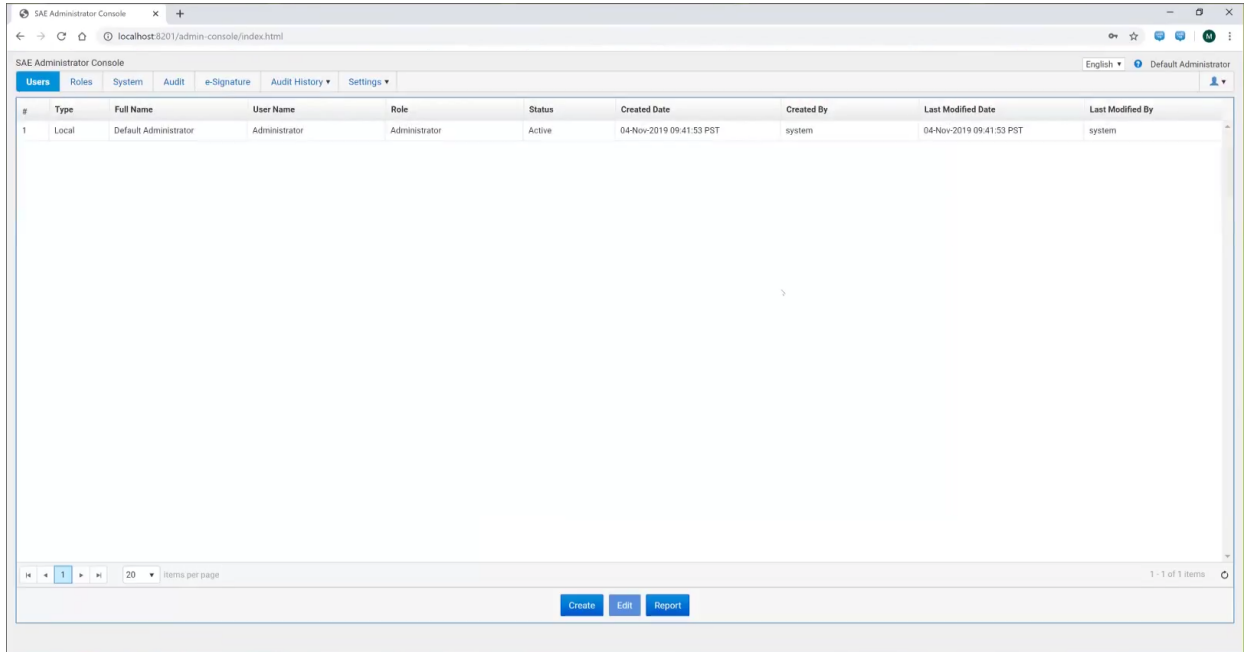
---

**Note:** If the SAE Administrator Console is configured to manage the **SAE module** for more than one application, you can create roles that specify permissions for more than one application (for example, Attune™ Cytometric Software and SeqStudio™ Genetic Analyzer). However, this section is focused solely on managing the **SAE module** for the Attune™ Cytometric Software.

---

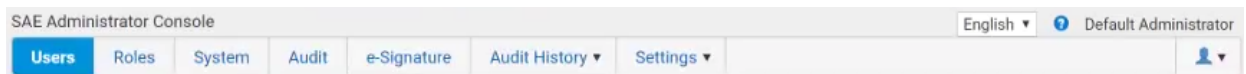
## Open the SAE Administrator Console

To open the SAE Administrator Console from the Attune™ Cytometric Software, click **View SAE Console** on the **SAE ribbon tab** (“Image Settings tab” on page 112). The SAE Administrator Console opens in the system’s default web browser.



## SAE Administrator Console tabs

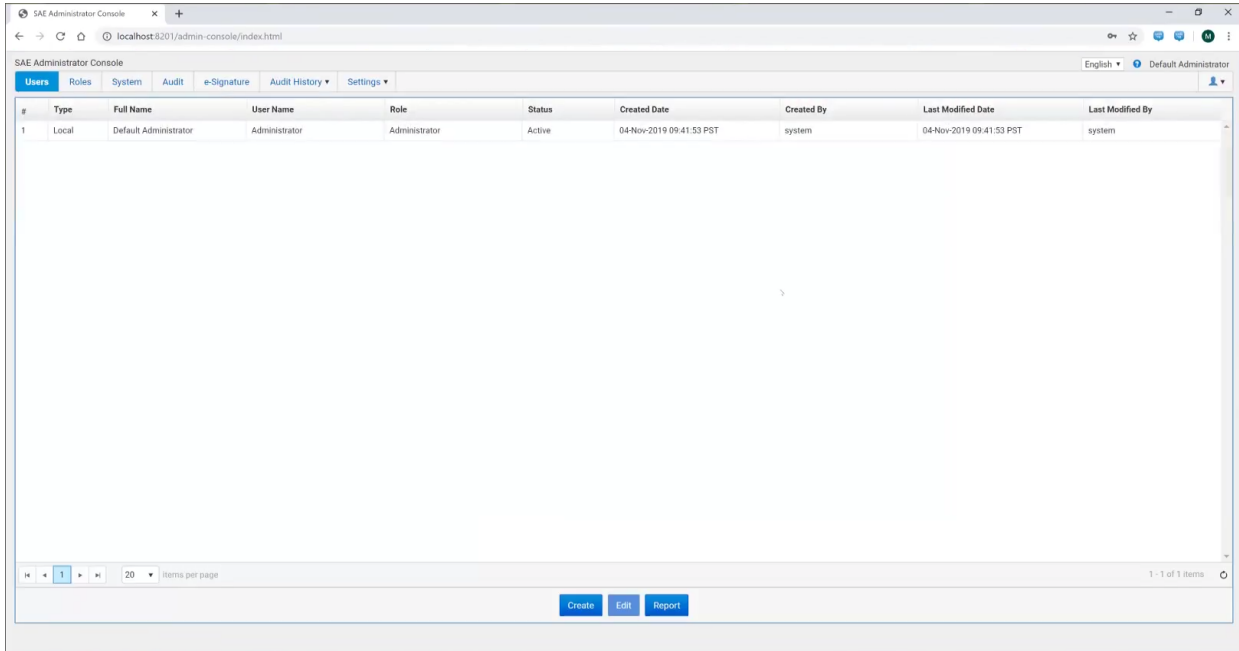
The SAE Administrator Console is organized into a series of tabs which represent the main functions of the **SAE Module**.



- **Users** (“Users tab” on page 882)
- **Roles** (“Roles tab” on page 886)
- **System** (“System tab” on page 897)
- **Audit** (“Audit tab” on page 903)
- **e-Signature** (“e-Signature tab” on page 909)
- **Audit History** (“Audit History tab” on page 913)
- **Settings** (“Settings tab” on page 927)

## Users tab

*Users tab* displays a list of the current SAE accounts in the SAE server, and allows the SAE Administrator to create additional SAE users, edit current users, and to create Users reports.



The screenshot shows the SAE Administrator Console interface. The 'Users' tab is selected, displaying a table with the following data:

#	Type	Full Name	User Name	Role	Status	Created Date	Created By	Last Modified Date	Last Modified By
1	Local	Default Administrator	Administrator	Administrator	Active	04-Nov-2019 09:41:53 PST	system	04-Nov-2019 09:41:53 PST	system

At the bottom of the table, there are three action buttons: **Create**, **Edit**, and **Report**. The page also shows a navigation bar with tabs for Users, Roles, System, Audit, e-Signature, Audit History, and Settings. The current user is identified as 'Default Administrator' in the top right corner.

- The list of SAE users contains columns for the User number (#), Type, Full Name, Username, Role, Status, Created Date, Created By, Last Modified Date, and Last Modified By.
- The full name of the currently sign in user is displayed on the upper right corner of the SAE Administrator Console.
- There are three action buttons in the Users tab:
  - **Create:** Allows the SAE Administrator to create additional SAE users (see “Create an SAE user account” on page 883).
  - **Edit:** Allows the SAE Administrator to edit the settings for the selected SAE user (see “Edit an SAE user account” on page 884).
  - **Report:** Allows the SAE Administrator to create a PDF report that lists the current SAE accounts saved in the SAE server (see “Print or view a user report” on page 885).



## Create an SAE user account

1. In the SAE Administrator Console main screen, click the **Users** tab.
2. Click **Create** to open the *Create User Account* dialog.

The screenshot shows the 'Create User Account' dialog box. It features a blue header with the text 'Create User Account'. Below the header, there are several input fields and dropdown menus. The fields are: 'User name', 'Password', 'Re-enter password', 'First name', 'MI', 'Last name', 'Phone', 'Email', and 'Comments'. There are also dropdown menus for 'Role' and 'Status' (set to 'Active'). A checkbox labeled 'User must set new password at next sign in' is checked. At the bottom right, there are 'Save' and 'Cancel' buttons.

3. Enter the **username, password, first name**, (optional) **middle initial**, and **last name**. The field limits are specified in the system security function settings (see “System tab” on page 897).

---

**Note:** First name, MI (middle initial), and last name are used to create the User Full Name, which is displayed as the name of the signed-in user. You cannot delete a user account or change the username after you save the user account.

---

4. (Optional) Deselect **User must set new password at next sign in**. By default, this option is selected. Users must specify a new password the first time they sign in to an application.

---

**Note:** The user account password automatically expires after the number of days that are specified in the system security function settings (see “System tab” on page 897).

---

5. Select the **Role** for the user account. Available default options are **Administrator, Advanced™ User, User, and Reviewer**.

The **No Privileges Role** is for internal use by the SAE Administrator Console. Do not assign this role to a user account.

---

**Note:** Each role grants specific SAE permissions to the user. You can also create a custom role with specific privileges in the Roles tab. For more information, see “Roles tab” on page 886.

---

6. Leave the **Status** set to **Active**.
7. (Optional) Enter phone, email (for information only), and comments.
8. Click **Save**. The newly created SAE user is displayed in the SAE users list in the Users tab.



## Edit an SAE user account

1. In the SAE Administrator Console main screen, click the **Users** tab.
2. Select a user account, then click **Edit** to open the *Edit User Account* dialog.
3. Edit the settings as desired. You cannot edit the username of an existing user.
4. Click **Save**.

---

**Note:** The Username cannot be changed after a user account is created and saved.

---

## Activate a suspended SAE user account

1. In the SAE Administrator Console main screen, click the **Users** tab.
2. Select a user account, then click **Edit** to open the *Edit User Account* window.
3. Change the Status from **Suspended** to **Active**.
4. Click **Save**.

## Disable (inactivate) an SAE user account

1. In the SAE Administrator Console main screen, click the **Users** tab.
2. Select a user account, then click **Edit** to open the *Edit User Account* window.
3. Change the Status from **Active** to **Inactive**.
4. Click **Save**.

## Reset an SAE user password

1. In the SAE Administrator Console main screen, click the **Users** tab.
2. Select a user account, then click **Edit** to open the *Edit User Account* window.
3. Enter a replacement password for the user account, then re-enter the password for confirmation.
4. If you assigned the user account a temporary password, then select **User must set new password at next sign in** to require the user to enter a new password at sign in.
5. Click **Save**.

---

**IMPORTANT!** There is no way to recover a forgotten password. If the SAE Administrator forgets their password, the software must be reinstalled. Export all data before reinstalling the software. Otherwise, the data will be lost. For more information, see “Export configuration” on page 932.

---



## Print or view a user report

1. In the SAE Administrator Console main screen, click the **Users** tab.
2. Click **Report**. The user report opens in the default web browser.
3. View the report, then print or save the report as a PDF in the desired location.
4. Close the report.

## Roles tab

**Roles** tab enables SAE Administrators to create, edit, and delete custom roles, and to create roles reports. SAE roles determine the SAE permissions that are associated with an SAE user account.

#	Role	Description	Users	Created Date	Created By	Last Modified Date	Last Modified By
1	Administrator	System role with full privileges	1	04-Nov-2019 09:41:52 PST	system	04-Nov-2019 09:41:52 PST	system
2	No Privileges Role	System role with no privileges	0	04-Nov-2019 09:41:52 PST	system	04-Nov-2019 09:41:52 PST	system
3	Advanced User	Attune - Advanced User	0	04-Nov-2019 09:41:57 PST	system	04-Nov-2019 09:41:57 PST	system
4	User	Attune - Basic User	0	04-Nov-2019 09:41:57 PST	system	04-Nov-2019 09:41:57 PST	system

- The **SAE module** for the Attune™ Cytometric Software has the following default SAE account types:
  - **Administrator:** System administrator role with full access to the software and full privileges, including the ability to modify SAE configuration and define the system security policy.
  - **Advanced User:** Similar to Attune™ advanced user.
  - **User:** Similar to Attune™ user.
  - **Reviewer:** Can only review **Audit History** and **Performance Test Reports**.
  - **No Privileges Role:** This role is for internal use only by the SAE Administrator Console to set up user repositories. Do not assign this role to a user account.
- For the permissions assigned to each default SAE account type, see “SAE account permissions” on page 887.
- There are four action buttons in the **Roles** tab:
  - **Create:** Enables the SAE Administrator to create custom SAE roles that give granular permissions to the Attune™ Cytometric Software features (see “Create a role” on page 894).
  - **Edit:** Enables the SAE Administrator to edit the settings for the selected custom SAE role (see “Edit a role” on page 895).  
The default SAE roles (**Administrator**, **Advanced User**, **User**, **Reviewer**, and **No Privileges**) are non-modifiable and are fixed with the functions they can perform.
  - **Delete:** Enables the SAE Administrator to delete selected custom SAE roles (see “Edit a role” on page 895).
  - **Report:** Enables the SAE Administrator to create a PDF report that lists the available SAE roles (see “Print or view a role report” on page 896).

**Note:** Changes made to a role in the SAE Administrator Console are reflected in the client application (i.e., Attune™ Cytometric Software) within 10 seconds.

## SAE account permissions

The permissions assigned to each default SAE account type are listed below. To create a custom role with specific privileges, see “Create a role” on page 894.

Permission	Administrator	Advanced User	User	Reviewer
<b>SAE Administrator Console permissions</b>				
<b>Security Configuration</b>				
- Configure Security and Auditing	Yes	No	No	No
<b>Audit History</b>				
- View Action Records	Yes	No	No	Yes
- View System Configuration Records	Yes	No	No	Yes
- View Application Object Records	Yes	No	No	Yes
- View Instrument Run Records	Yes	No	No	Yes
<b>Attune™ Cytometric Software permissions</b>				
<b>Acquisition Control</b>				
- Copy and paste run protocols settings	Yes	Yes	Yes	No
- Modify collect options	Yes	Yes	Yes	No
- Modify run protocol settings	Yes	Yes	Yes	No
- Record over (overwrite) existing sample data	Yes	Yes	Yes	No
- Run and record samples/plates	Yes	Yes	Yes	No
<b>Compensation</b>				
- Apply compensation	Yes	Yes	Yes	No
- Create compensation	Yes	Yes	Yes	No
- Delete compensation	Yes	Yes	Yes	No
- Edit compensation channels	Yes	Yes	Yes	No
- Export compensation	Yes	Yes	Yes	No
- Import compensation	Yes	Yes	Yes	No
- Modify compensation values	Yes	Yes	Yes	No
<b>Experiment Management</b>				

(continued)

Permission	Administrator	Advanced User	User	Reviewer
- Change experiment and group colors	Yes	Yes	Yes	No
- Create experiments	Yes	Yes	Yes	No
- Create experiments from templates	Yes	Yes	Yes	No
- Create samples and groups	Yes	Yes	Yes	No
- Create templates	Yes	Yes	Yes	No
- Delete experiments	Yes	Yes	Yes	No
- Delete samples	Yes	Yes	Yes	No
- Delete templates	Yes	Yes	Yes	No
- Duplicate experiments	Yes	Yes	Yes	No
- Export experiments	Yes	Yes	Yes	No
- Export FCS/ACS files	Yes	Yes	Yes	No
- Export images	Yes	Yes	Yes	No
- Export sample list	Yes	Yes	Yes	No
- Export templates	Yes	Yes	Yes	No
- Import experiments	Yes	Yes	Yes	No
- Import FCS/ACS files	Yes	Yes	Yes	No
- Import sample list	Yes	Yes	Yes	No
- Import templates	Yes	Yes	Yes	No
- Modify experiment annotation	Yes	Yes	Yes	No
- Modify templates	Yes	Yes	Yes	No
- Open experiments	Yes	Yes	Yes	No
- Remove FCS files	Yes	Yes	Yes	No
- Show/Hide experiment and group colors	Yes	Yes	Yes	No
- Update keywords in exported FCS files	Yes	Yes	Yes	No
<b>Filter Configuration</b>				
- Create and edit filters	Yes	Yes	No	No
- Manage filter configurations	Yes	Yes	No	No
- Modify filter mapping	Yes	Yes	No	No



(continued)

Permission	Administrator	Advanced User	User	Reviewer
- View filter configuration	Yes	Yes	Yes	No
<b>Image Analysis Settings</b>				
- Allow Image Adjustment	Yes	Yes	Yes	No
- Apply Image Filtering	Yes	Yes	Yes	No
- Delete Cell Images	Yes	Yes	Yes	No
- Export Cell Images	Yes	Yes	Yes	No
- Insert Cell Image Objects	Yes	Yes	Yes	No
- Measure Image	Yes	Yes	Yes	No
- Modify Mask Settings	Yes	Yes	Yes	No
- Show/Hide Masks	Yes	Yes	Yes	No
<b>Image Capture Settings</b>				
- Modify camera control settings	Yes	Yes	Yes	No
- Modify camera ROI settings	Yes	Yes	Yes	No
- Modify image capture settings	Yes	Yes	Yes	No
<b>Image Processing</b>				
- Overwrite image processing	Yes	Yes	Yes	No
- Process images	Yes	Yes	Yes	No
- Remove image processing data	Yes	Yes	Yes	No
<b>Instrument Control</b>				
- Run auto sampler calibration	Yes	Yes	Yes	No
- Run debubble and unclog	Yes	Yes	Yes	No
- Run deep clean and shutdown	Yes	Yes	Yes	No
- Run recover sample	Yes	Yes	Yes	No
- Run rinse, SIP sanitize	Yes	Yes	Yes	No
- Run startup	Yes	Yes	Yes	No
- Run system decontamination	Yes	Yes	No	No
- Run system self test	Yes	Yes	No	No
<b>Instrument Settings</b>				

(continued)

Permission	Administrator	Advanced User	User	Reviewer
- Copy and paste instrument settings	Yes	Yes	Yes	No
- Export instrument settings	Yes	Yes	Yes	No
- Import instrument settings	Yes	Yes	Yes	No
- Modify advanced instrument settings	Yes	Yes	No	No
- Modify instrument configuration	Yes	Yes	No	No
- Modify parameter on/off states	Yes	Yes	Yes	No
- Modify parameter target and label names	Yes	Yes	Yes	No
- Modify system instrument settings	Yes	No	No	No
- Modify thresholds	Yes	Yes	Yes	No
- Modify voltages	Yes	Yes	Yes	No
<b>Options</b>				
- Create and edit plates	Yes	Yes	No	No
- Create and edit user keywords	Yes	Yes	Yes	No
- Manage global keywords	Yes	No	No	No
- Modify administrator options	Yes	No	No	No
- Modify configuration options	Yes	No	No	No
- Modify default colors	Yes	Yes	Yes	No
- Modify default font options	Yes	Yes	Yes	No
- Modify default gate options	Yes	Yes	Yes	No
- Modify default plot options	Yes	Yes	Yes	No
- Modify default sample and group name	Yes	Yes	Yes	No
- Modify default statistics options	Yes	Yes	Yes	No
- Modify display options	Yes	Yes	Yes	No
- Modify export options	Yes	Yes	Yes	No
- Modify images options	Yes	No	No	No
<b>Performance Test</b>				
- Reset baseline	Yes	Yes	No	No
- Run performance test	Yes	Yes	No	No





(continued)

Permission	Administrator	Advanced User	User	Reviewer
- View performance test report	Yes	Yes	Yes	Yes
<b>Workspace and Overlay Gates</b>				
- Change gate color	Yes	Yes	Yes	No
- Change gate coordinates	Yes	Yes	Yes	No
- Change gate name	Yes	Yes	Yes	No
- Change gate opacity	Yes	Yes	Yes	No
- Change gate type	Yes	Yes	Yes	No
- Create gates	Yes	Yes	Yes	No
- Delete gates	Yes	Yes	Yes	No
- Export gate to FCS file	Yes	Yes	Yes	No
- Modify autogate settings	Yes	Yes	Yes	No
- Modify gate equation	Yes	Yes	Yes	No
- Modify gate Z order	Yes	Yes	Yes	No
- Show/Hide gate name	Yes	Yes	Yes	No
<b>Workspace and Overlay Plots</b>				
- Change density plot color and mode	Yes	Yes	Yes	No
- Change legend text	Yes	Yes	Yes	No
- Change overlay opacity	Yes	Yes	Yes	No
- Change overlay plot color	Yes	Yes	Yes	No
- Change percent of displayed events	Yes	Yes	Yes	No
- Change plot axes labels	Yes	Yes	Yes	No
- Change plot parameters	Yes	Yes	Yes	No
- Change plot range	Yes	Yes	Yes	No
- Change plot resolution	Yes	Yes	Yes	No
- Change plot scale	Yes	Yes	Yes	No
- Change plot title	Yes	Yes	Yes	No
- Change plot types	Yes	Yes	Yes	No
- Create overlays	Yes	Yes	Yes	No

(continued)

Permission	Administrator	Advanced User	User	Reviewer
- Delete overlays	Yes	Yes	Yes	No
- Delete plots	Yes	Yes	Yes	No
- Format plot text	Yes	Yes	Yes	No
- Hide tick marks on overlays	Yes	Yes	Yes	No
- Insert plots	Yes	Yes	Yes	No
- Modify histogram properties	Yes	Yes	Yes	No
- Modify overlay 3D options	Yes	Yes	Yes	No
- Modify plot statistics	Yes	Yes	Yes	No
- Perform overlay calculations	Yes	Yes	Yes	No
- Reorder overlay plot members	Yes	Yes	Yes	No
- Save plot as image	Yes	Yes	Yes	No
- Show/Hide overlay plots	Yes	Yes	Yes	No
- Show legend on overlays	Yes	Yes	Yes	No
<b>Workspace Layout</b>				
- Align objects	Yes	Yes	Yes	No
- Change workspace grid size	Yes	Yes	Yes	No
- Modify workspace mode	Yes	Yes	Yes	No
- Move and resize objects	Yes	Yes	Yes	No
<b>Other Workspace Objects</b>				
- Copy objects to clipboard	Yes	Yes	Yes	No
- Delete images	Yes	Yes	Yes	No
- Delete statistics	Yes	Yes	Yes	No
- Delete text boxes	Yes	Yes	Yes	No
- Edit text box text	Yes	Yes	Yes	No
- Format statistics	Yes	Yes	Yes	No
- Format text box	Yes	Yes	Yes	No
- Insert images	Yes	Yes	Yes	No
- Insert statistics	Yes	Yes	Yes	No



(continued)

Permission	Administrator	Advanced User	User	Reviewer
- Insert text boxes	Yes	Yes	Yes	No
- Modify statistics box statistics	Yes	Yes	Yes	No
<b>Other</b>				
- Export workspace	Yes	Yes	Yes	No
- Import workspace	Yes	Yes	Yes	No
- Modify heatmap settings	Yes	Yes	Yes	No
- Printing and page layout	Yes	Yes	Yes	No
- View system logs	Yes	Yes	Yes	No
- View user logs	Yes	No	No	No
<b>Security Configuration</b>				
- Perform e-signing	Yes	Yes	No	No

## Create a role

1. In the SAE Administrator Console main screen, click the **Roles** tab.
2. Click **Create** to open the **Create Role dialog**.

**Create Role**

\* Name

Description

**Permissions**

- SAE Administrator Console
  - Security Configuration
  - Audit History
- Attune Software
  - Acquisition Control
  - Compensation
  - Experiment Management
  - Filter Configuration
  - Image Analysis Settings
  - Image Capture Settings
  - Instrument Control
  - Instrument Settings
  - Options
  - Performance Test
  - Workspace and Overlay Gates
  - Workspace and Overlay Plots
  - Workspace Layout
  - Other Workspace Objects
  - Other
  - Security Configuration

3. Enter the **Name** and (*optional*) **Description** for the new role.
4. Select the SAE **Permissions** for the role.  
To select all **Permissions** in a category, select the **Permissions checkbox** next to the category.
5. Click **Save**. The newly created role is displayed in the **SAE Roles list** in the **Roles** tab and becomes available in the **Role dropdown** in the **Create User Account dialog** (“Create an SAE user account” on page 883).

## Edit a role

1. In the SAE Administrator Console main screen, click the **Roles** tab.
2. Select a role, then click **Edit** to open the **Edit Role dialog**.

The screenshot shows the 'Edit Role' dialog box. The 'Name' field is set to 'TestRole'. The 'Description' field is empty. The 'Permissions' section is expanded, showing a tree view of permissions. Under 'SAE Administrator Console', 'Security Configuration' and 'Audit History' are unchecked. Under 'Attune Software', several permissions are checked, including 'Acquisition Control', 'Compensation', 'Experiment Management', 'Filter Configuration', 'Image Analysis Settings', 'Image Capture Settings', 'Instrument Control', 'Instrument Settings', 'Options', 'Performance Test', 'Workspace and Overlay Gates', 'Workspace and Overlay Plots', 'Workspace Layout', 'Other Workspace Objects', 'Other', and 'Security Configuration'. At the bottom right, there are 'Save' and 'Cancel' buttons.

---

**Note:** You cannot edit the default SAE roles (**Administrator**, **Advanced User**, **User**, **Reviewer**, and **No Privileges**).

---

3. Edit the settings as desired, then click **Save**.

---

**Note:** Changes made to a role in the SAE Administrator Console are reflected in the client application (i.e., Attune™ Cytometric Software) within 10 seconds.

---

## Delete a role

1. In the SAE Administrator Console main screen, click the **Roles** tab.
2. Select a role, then click **Delete**.

---

**Note:** If an SAE user account is assigned to a role, that role cannot be deleted.

---



## Print or view a role report

1. In the SAE Administrator Console main screen, click the **Roles** tab.
2. Click **Report**. The role report opens in the default web browser.
3. View the report, then print or save the report as a PDF in the desired location.
4. Close the report.

## System tab

**System** tab enables the **SAE Administrator** to modify SAE configuration and define the system security policy. In this tab, the SAE Administrator can:

- Configure username settings (“User Name Settings” on page 897)
- Set password policy (“Password Policy” on page 898)
- Define account lockout policy (“Account Lockout Policy” on page 899)
- Configure other system settings (“Other Settings” on page 900).

Access to the **System** tab is restricted to SAE users with the permission to **Configure Security and Auditing** (see “SAE account permissions” on page 887).

The screenshot shows the SAE Administrator Console interface. The top navigation bar includes 'Users', 'Roles', 'System' (selected), 'Audit', 'e-Signature', 'Audit History', and 'Settings'. The main content area is titled 'Password Policy' and contains the following settings:

- Minimum length: 8 characters
- Maximum length: 32 characters
- May not reuse previous: 3 passwords
- Password complexity: Alphabets (Allowed), Uppercase (Allowed), Lowercase (Allowed), Numeric (Allowed), Special characters (Allowed)
- Maximum password age: 120 days
- Minimum password age: 0 days
- Password expiry reminder: Enabled
- Send reminder: 3 days
- Username check: Enabled
- Compromised phrase check: Enabled

At the bottom of the form, there are two buttons: 'Reset to Defaults' and 'Apply Settings'.

**Note:** Settings in the **System** tab affect all SAE user accounts. Settings are applied the next time that users sign in to an application.

## User Name Settings

**User Name Settings** enable the SAE Administrator to set a minimum and maximum length for SAE user names.

The screenshot shows the 'User Name Settings' form with the following settings:

- Minimum length: 8 characters
- Maximum length: 32 characters

- By default, minimum and maximum length for SAE user names are set as 8 and 32 characters, respectively.
- The Attune™ Cytometric Software only allows a maximum of 50 characters for user names. If a user name that is longer than 50 characters is chosen, the SAE user cannot login into the Attune™ Cytometric Software.

## Password Policy

**Password Policy** provides options to set up minimum and maximum password length, password complexity requirements (inclusion of alpha, numeric, lower case, upper case and special characters), password age, password expiration reminder, username check, and compromised phrase check settings.

Password Policy

<b>Minimum length</b> <input style="width: 60px;" type="text" value="8"/> characters	<b>Maximum length</b> <input style="width: 60px;" type="text" value="64"/> characters
<b>May not reuse previous</b> <input style="width: 60px;" type="text" value="3"/> passwords	
<b>Password complexity</b>	
<b>Alphabets</b> <input type="radio"/> Forbidden <input checked="" type="radio"/> Allowed	<b>at least</b> <input style="width: 40px;" type="text" value="0"/> occurrences
<b>Uppercase</b> <input type="radio"/> Forbidden <input checked="" type="radio"/> Allowed	<b>at least</b> <input style="width: 40px;" type="text" value="0"/> occurrences
<b>Lowercase</b> <input type="radio"/> Forbidden <input checked="" type="radio"/> Allowed	<b>at least</b> <input style="width: 40px;" type="text" value="0"/> occurrences
<b>Numeric</b> <input type="radio"/> Forbidden <input checked="" type="radio"/> Allowed	<b>at least</b> <input style="width: 40px;" type="text" value="0"/> occurrences
<b>Special characters</b> <input type="radio"/> Forbidden <input checked="" type="radio"/> Allowed	<b>at least</b> <input style="width: 40px;" type="text" value="0"/> occurrences
<b>Maximum password age</b> <input style="width: 60px;" type="text" value="120"/> days	<b>Minimum password age</b> <input style="width: 60px;" type="text" value="0"/> days
<b>Password expiry reminder</b> <input checked="" type="radio"/> Enabled <input type="radio"/> Disabled	<b>Send reminder</b> <input style="width: 60px;" type="text" value="3"/> days
<b>Username check</b> <input checked="" type="radio"/> Enabled <input type="radio"/> Disabled	<b>Compromised phrase check</b> <input checked="" type="radio"/> Enabled <input type="radio"/> Disabled

- The password policy is only applied to SAE user accounts and is considered when a user chooses to change their password in the Attune™ Cytometric Software (see “Change a password” on page 712) or is required to change their password on login (see “Expired password” on page 41).
- If the password does not meet any of the password policy settings, a balloon tip is displayed specifying the password policy criteria.
- **Password Age** settings specify the maximum (password expiration) and minimum password age (time before a user can change their password), and the password expiration reminder settings.
- When the password reminder setting is active, the user is presented with a warning dialog when logging in telling them that their password will expire in a specified number of days (as set by the send reminder value).



- **Username check**, when selected, searches for user identifiable and easy to guess phrases in the chosen password. If such phrases are found, the software displays a prompt for the user to choose a password that is not based on the username and that does not contain multiple sequential numbers.
- **Compromised phrase check**, when selected, searches for commonly used word or phrase in the chosen password. If such a word or phrase is found in the password, the software displays a prompt for the user to select a different password.

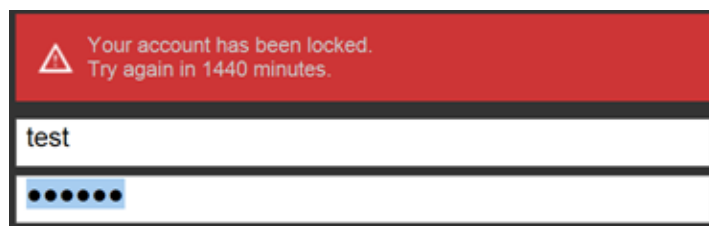
## Account Lockout Policy

**Account Lockout Policy** settings control the number of failed sign in attempts allowed as an SAE user before locking the user out for a specified time.

The screenshot shows the 'Account Lockout' configuration page. At the top, there is a header 'Account Lockout'. Below it, there are four main settings:

- Account lockout:** A radio button group with 'Enabled' selected and 'Disabled' unselected.
- Threshold:** A text input field containing the number '3', with a 'sign in attempts' label to its right.
- Account lockout duration:** A text input field containing '1440', with a 'minutes' label to its right.
- Sign in attempts counter reset:** A radio button group with 'Enabled' selected and 'Disabled' unselected.
- Reset failure sign in counter after:** A text input field containing '1440', with a 'minutes' label to its right.

- By default, the **Account Lockout** is enabled and allows 3 invalid sign in attempts before locking the account (suspending) for a specified time (24 hours by default). You can set the lockout threshold from 1 to 999 attempts, and the lockout time from 1 to 99,999 minutes.
- The account lockout policy also includes a setting that allows the sign in attempt counter to automatically reset (enabled by default) after a specified time (24 hours by default and configurable between 1 and 99,999 minutes).
- If a user exceeds the sign in attempt limit, the user's account is suspended and a warning is displayed in the sign in screen.



## Other Settings

**Other Settings** enables the **SAE Administrator** to control the following settings:

- Automatic screen locking (due to inactivity) (“Other Settings” on page 900)
- Open files from non-SAE system (“Open file from non-SAE system” on page 901)
- Client offline sign in (“Client offline sign in” on page 901)



Other Settings

Automatic screen locking  Enabled  Disabled

Inactivity duration 30 minutes

Open file from non-SAE system  Forbidden  Allowed

Client offline sign in  Enabled  Disabled

Offline sign in threshold 0 minutes

### Automatic screen locking

**Automatic screen locking** allows the system to automatically lock out users after a specified period of inactivity, which is defined as no keyboard or mouse input in the Attune™ Cytometric Software.

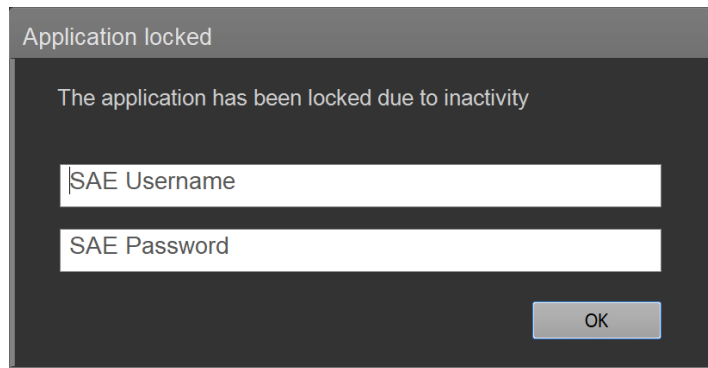
- Automatic screen locking is enabled by default and set to 30 minutes of inactivity. You can set the inactivity length from 1 to 99,999 minutes.
- When you sign in as an SAE user, your session is actively managed to ensure that the system is in use by you. The software registers any movement of the mouse or keyboard input as activity.

---

**Note:** You can also manually lock the Attune™ Cytometric Software by selecting **Lock Application** from the **File** menu or by entering **Ctrl+Alt+L** on the keyboard.

---

- When the SAE server is configured to automatically lock out users after a specified period of inactivity, the **Application Locked dialog** is displayed after the set period of inactivity, requiring you to enter credentials to unlock the software.



- During acquisition, when the lockout dialog is displayed due to inactivity, the acquisition continues uninterrupted.
- If you are signed in when the server is offline, automatic account lockout does not take effect.

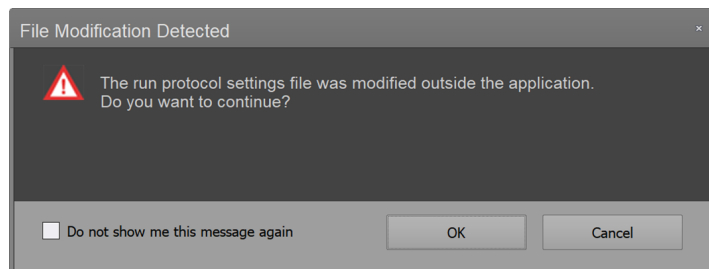
## Open file from non-SAE system

When allowed, **Open file from non-SAE system** enables you to access and open files (such as FCS files, Workspace files, Run protocol files) from non-SAE systems. This feature also enables you to open, duplicate, and export Experiments containing modifications that result in audit gaps.

**Note:** When an Experiment is opened, all associated files (such as FCS files, Workspace files, Run protocol files) are checked for any modifications by calculating file checksums and comparing them to the expected values in the database. A mismatch between the checksums from the opened file and the expected values results in an audit gap. When an audit gap is detected, the software displays the **File Modification Detected** warning dialog (see below).

- By default, **Open file from non-SAE system** is set to **Forbidden**.
- To transfer data from a non-SAE Attune™ account into an SAE account, set **Open file from non-SAE system** to **Allowed**.
- For a first migration step for files that do not have the SAE marking, the **SAE Administrator** can allow non-SAE files to be opened for set time period, then readjust the SAE marking requirement to not allow those files after the first migration.
- When set to **Allowed**, **Open file from non-SAE system** also enables you to open an Experiment that has been altered outside of the application and continue with that Experiment. However, depending on what has been changed in the Experiment data file, continuing with a tampered Experiment can result in an Audit gap.

If there are audit gaps in the audit history, the Attune™ Cytometric Software displays the **File Modification Detected** warning dialog to notify you that you are trying to open an Experiment that has an invalid audit history and ask you to confirm to continue.

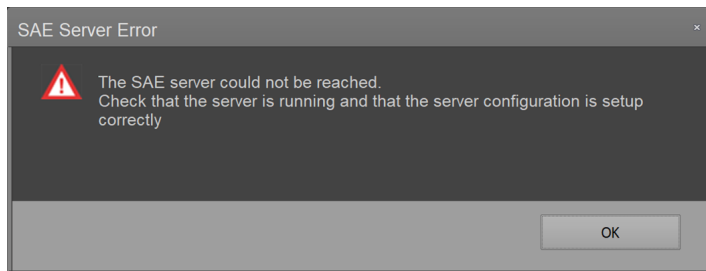


## Client offline sign in

**Client offline sign in** settings control whether an SAE user can sign in the Attune™ Cytometric Software when the SAE server is offline or disconnected.

- **Client offline sign in** is only available for SAE accounts that have signed into the software at least once when the server was online.  
If an account has not accessed the Attune™ Cytometric Software before, that account cannot perform offline sign in, even if the **Client offline sign in** option is enabled for the system.
- By default, **Client offline sign in** is disabled and set to allow an offline login threshold of 0 minutes.
- The **Offline sign in threshold** specifies the required elapsed time since the last disconnected state with the SAE server and when an SAE user can sign in the Attune™ Cytometric Software.  
When **Client offline sign in** is enabled, a value between 1 and 99,999 minutes must be specified for the Offline sign in threshold.

- If the **Offline sign in threshold** is exceeded, offline sign in cannot be performed, and the software displays the **SAE Server Error** message.

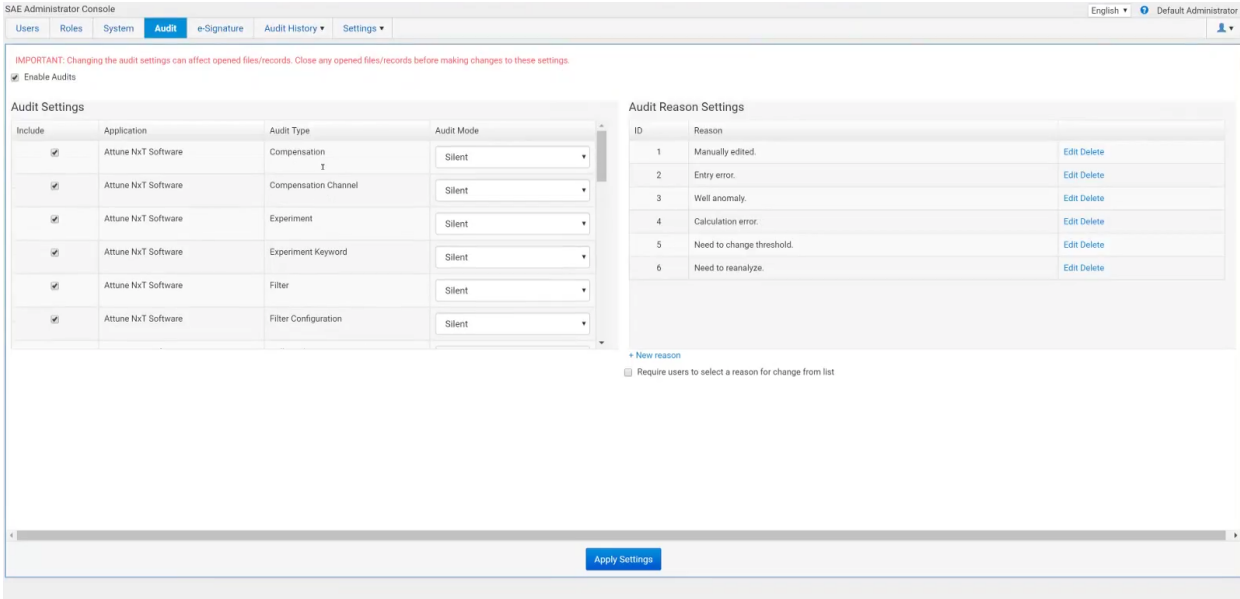


- A currently signed in user can remain signed in after the sign in threshold limit is exceeded.
- After an SAE user has signed in offline, the security configuration (FAC) that takes effect is based on the last connected state.
- Audited events, data audits and e-Signature records are generated and sent to server when the SAE server connection is restored. The audit data and e-Signature data are stored in an encrypted cache and synchronized with the server on reconnection.
- Automatic account lockout, password reminder notification, and mandatory password change does not take effect under offline sign in.
- The following functions cannot be performed under **Client offline sign in**:
  - **Disable SAE**: If a user attempts to disable the SAE mode when in the offline mode, the software displays the **SAE Server Error** message stating that the SAE mode cannot be disabled when in the offline mode.
  - **Change Password (for SAE user)**: If a user attempts to change the password when in the offline mode, the software displays the **SAE Server Error** message stating that the password cannot be changed when in the offline mode.

## Audit tab

**Audit** tab allows you to enable or disable audits, which track actions performed by users, changes to SAE settings, and changes made to application objects (for example, Experiments). In the **Audit** tab, you can also control which actions and application objects are audited, and require users to provide a reason for changes made to auditable application objects.

**IMPORTANT!** Changing audit settings after Experiments have been created can result in audit gaps. Ensure that the audit settings are configured before using the Attune™ Cytometric Software in the SAE mode.



SAE Administrator Console

English • Default Administrator

Users Roles System **Audit** e-Signature Audit History Settings

**IMPORTANT:** Changing the audit settings can affect opened files/records. Close any opened files/records before making changes to these settings.

Enable Audits

**Audit Settings**

Include	Application	Audit Type	Audit Mode
<input checked="" type="checkbox"/>	Attune NXT Software	Compensation	Silent
<input checked="" type="checkbox"/>	Attune NXT Software	Compensation Channel	Silent
<input checked="" type="checkbox"/>	Attune NXT Software	Experiment	Silent
<input checked="" type="checkbox"/>	Attune NXT Software	Experiment Keyword	Silent
<input checked="" type="checkbox"/>	Attune NXT Software	Filter	Silent
<input checked="" type="checkbox"/>	Attune NXT Software	Filter Configuration	Silent

**Audit Reason Settings**

ID	Reason	
1	Manually edited.	Edit Delete
2	Entry error.	Edit Delete
3	Well anomaly.	Edit Delete
4	Calculation error.	Edit Delete
5	Need to change threshold.	Edit Delete
6	Need to reanalyze.	Edit Delete

+ New reason

Require users to select a reason for change from list

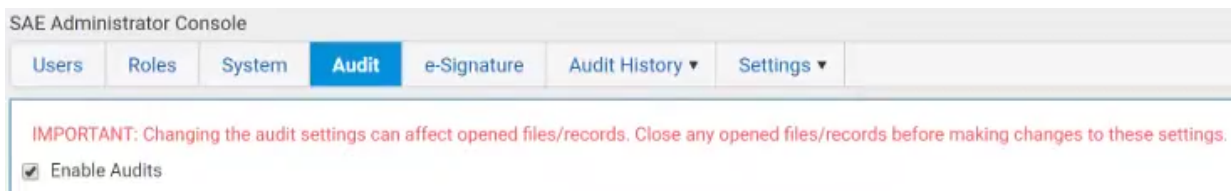
Apply Settings

**Audit** tab has the following two functional groups:

- **Audit Settings:** Consists of the **Enable Audits** option (“Enable Audits” on page 904) and the **Audit Settings** panel (“Audit Settings panel” on page 905)
- **Audit Reason Settings** (“Audit Reason Settings panel” on page 906)

## Enable Audits

**Enable Audits** option turns the audit function on or off (see “Enable or disable the audit function” on page 907).



- When the SAE mode is active and **Enable Audits** option in the **Audit** tab is selected, any change to an Experiment results in an audit transaction that is captured in the SAE console, where the old and new values of each Experiment attribute is recorded.
- You can disable the audit function in its entirety or for specific **Audit Types** (see “Audit Settings panel” on page 905).
- By default, auditing is active for all **Audit Types**.
- If you have permission to view **Audit History**, you can view the audit records in the **Audit History** tab (“Audit History tab” on page 913).

---

**Note:** You cannot disable auditing for **Action Records** (“Action Records audit log” on page 914) and **System Configuration** (“System Configuration audit log” on page 917). Even if you unselect the **Enable Audits** checkbox, these logs are always audited.

---



## Audit Settings panel

**Audit Settings** panel enables you to select or deselect which specific application objects are included in the audit, and the **Audit mode** (see “Select items to audit and set the audit mode” on page 907).

Include	Application	Audit Type	Audit Mode
<input checked="" type="checkbox"/>	Attune NxT Software	Compensation	Silent
<input checked="" type="checkbox"/>	Attune NxT Software	Compensation Channel	Silent
<input checked="" type="checkbox"/>	Attune NxT Software	Experiment	Silent
<input checked="" type="checkbox"/>	Attune NxT Software	Experiment Keyword	Silent
<input checked="" type="checkbox"/>	Attune NxT Software	Filter	Silent
<input checked="" type="checkbox"/>	Attune NxT Software	Filter Configuration	Silent

- **Audit Settings** panel has four columns:
  - **Include:** Selects the items to include in the audit.
  - **Application:** Displays the application to which the audit applies.  
If the **SAE Administrator Console** is configured to manage the SAE module for more than one application, you can audit more than one application (for example, Attune™ Cytometric Software and SeqStudio™ Genetic Analyzer).
  - **Audit Type:** Lists what is included in the audit.
  - **Audit Mode:** Enables you to select one of three audit modes:

Audit Mode	Reason
<b>Silent</b>	The event is audited, no reason prompt is displayed.
<b>Optional</b>	The event is audited, a reason prompt is displayed, but you can cancel and continue without entering a reason.
<b>Required</b>	The event is audited, a reason prompt is displayed, and you must specify a reason.

- The default **Audit Mode** for Attune™ Cytometric Software application objects is **Silent**.
- In **Silent** auditing, audits take place silently with no user input. The changes are audited with the following default reasons (depending on **Change** type):

Change type	Reason
Create	User wants to create
Update	User wants to update
Delete	User wants to delete

- If an application object is set to audit using the **Optional** or **Required** mode, the **Enter Audit Reason** dialog is displayed each time a change is made that results in an audit.

## Audit Reason Settings panel

**Audit Reason Settings** panel enables you to configure the audit reasons that are presented in the **Enter Audit Reason** dialog.

### Audit Reason Settings

ID	Reason	
1	Manually edited.	<a href="#">Edit</a> <a href="#">Delete</a>
2	Entry error.	<a href="#">Edit</a> <a href="#">Delete</a>
3	Well anomaly.	<a href="#">Edit</a> <a href="#">Delete</a>
4	Calculation error.	<a href="#">Edit</a> <a href="#">Delete</a>
5	Need to change threshold.	<a href="#">Edit</a> <a href="#">Delete</a>
6	Need to reanalyze.	<a href="#">Edit</a> <a href="#">Delete</a>

[+ New reason](#)

- The SAE console is preloaded with the following defined audit reasons:
  - Manually edited
  - Entry error
  - Well anomaly
  - Calculation error
  - Need to change threshold
  - Need to reanalyze
- You can edit or delete any existing reason by clicking **Edit** or **Delete**.
- You can create new audit reasons by clicking **New reason**.



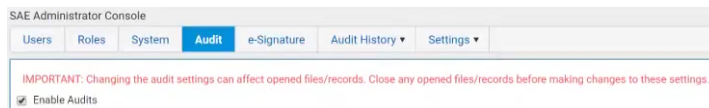


- You can enforce the use of predefined reasons instead of allowing a user to manually type a reason by selecting the **Require users to select a reason for change from list** option.

**Require users to select a reason for change from list**

## Enable or disable the audit function

- In the SAE Administrator Console main screen, click the **Audit** tab.
- Select or deselect **Enable Audits**. The role report opens in the default web browser.




---

**IMPORTANT!** Changing the audit settings can affect opened files or records. Close any opened files or records before making changes to audit settings.

---

- (Optional)* Set or modify the **Audit Settings** and the **Audit Reason Settings**.

---

**Note:** When the **Audit mode** is set to **Silent**, audit reasons are not available for user selection in an application.

---

- Click **Apply Settings**.

---

**Note:** You cannot turn off auditing for Action Records (“Action Records audit log” on page 914) and System Configuration (“System Configuration audit log” on page 917). Even if you unselect the **Enable Audits** checkbox, these are always audited.

---

## Select items to audit and set the audit mode

- In the SAE Administrator Console main screen, click the **Audit** tab.
- In the **Audit Settings** panel, select the items to audit.

Include	Application	Audit Type	Audit Mode
<input checked="" type="checkbox"/>	Attune NxT Software	Compensation	Silent
<input checked="" type="checkbox"/>	Attune NxT Software	Compensation Channel	Silent
<input checked="" type="checkbox"/>	Attune NxT Software	Experiment	Silent
<input checked="" type="checkbox"/>	Attune NxT Software	Experiment Keyword	Silent
<input checked="" type="checkbox"/>	Attune NxT Software	Filter	Silent
<input checked="" type="checkbox"/>	Attune NxT Software	Filter Configuration	Silent

3. Select the **Audit Mode** for each item you include for auditing:
  - **Silent:** The event is audited, no reason prompt is displayed.
  - **Optional:** The event is audited, a reason prompt is displayed, but the user can cancel and continue without entering a reason.
  - **Required:** The event is audited, a reason prompt is displayed, and the user must specify a reason.
4. Click **Apply Settings**.

## Configure audit reason settings

1. In the SAE Administrator Console main screen, click the **Audit** tab.  
Configure the audit reason settings in the Audit Reason Settings panel as described below.

Audit Reason Settings

ID	Reason	
1	Manually edited.	<a href="#">Edit</a> <a href="#">Delete</a>
2	Entry error.	<a href="#">Edit</a> <a href="#">Delete</a>
3	Well anomaly.	<a href="#">Edit</a> <a href="#">Delete</a>
4	Calculation error.	<a href="#">Edit</a> <a href="#">Delete</a>
5	Need to change threshold.	<a href="#">Edit</a> <a href="#">Delete</a>
6	Need to reanalyze.	<a href="#">Edit</a> <a href="#">Delete</a>

[+ New reason](#)

Require users to select a reason for change from list

2. (Optional) Select **Require users to select a reason for change from list** to require users to select a pre-defined audit reason from the Reason list.

### Add new audit reason

1. Click **New Reason**.
2. Enter a reason for change, then click **Save**.
3. Click **Apply Settings**.

### Edit an existing audit reason:

1. Click **Edit**.
2. Edit the reason for change, then click **Save**.
3. Click **Apply Settings**.

### Delete an existing audit reason:

1. Click **Delete**.
2. Click **Delete** to confirm deletion of the audit reason or **Cancel** to exit the dialog.
3. Click **Apply Settings**.

---

**Note:** Deleting an audit reason also deletes its ID number. The deleted ID number is not reused for the next audit reason in the list.

---



## e-Signature tab

**e-Signature** tab enables you to control the e-Signature rights of SAE roles, the actions requiring signatures, the reasons available for e-Signatures, and the data to be signed.

SAE Administrator Console

English | Default Administrator

Users | Roles | System | Audit | **e-Signature** | Audit History | Settings

**IMPORTANT:** Changing the e-Signature settings can affect opened files/records. Close any opened files/records before making changes to these settings.

Enable e-Signatures

Show e-signature configuration for: Attune NxT Software

**e-Signature Meanings**

Meanings

- Review and Approve Experiment for Template
- Review and Approve Data

+ New meaning

**Data signed for selected meaning**

Sign:  Data: Tube or Plate Experiment

**Actions Requiring Signatures**

Include	Action
<input type="checkbox"/>	Export Data
<input type="checkbox"/>	Print Data
<input type="checkbox"/>	Save as Template

**Number of signatures required for selected action**

Meanings	Administrator	Advanced User	User
Review and Approve Experiment for Template	0	0	0
Review and Approve Data	0	0	0

Apply Settings

- Actions that require signatures to be completed cannot be performed unless the required signatures are provided.
- Signatures represent the state of the Experiment at the time of signing. If any modifications are made to the Experiment after signing, actions that require signatures cannot be performed unless new signatures are provided.
- You can view the log of e-Signatures and the status of the signatures (Current or Obsolete) in the **Audit History** ▶ **Application Object Records** ▶ **e-Signature Records** tab (“e-Signature Records tab” on page 923).
- You can enable or disable the requirement for e-Signatures by selecting or unselecting the **Enable e-Signatures** checkbox (see “Enable or disable the e-Signature function” on page 910).
- There are three built-in **Actions Requiring Signatures** for the Attune™ Cytometric Software that you can select in **e-Signature** tab:
  - **Export Data**
  - **Print Data**
  - **Save As Template**

By default, the **Include** checkboxes for these actions are unchecked (that is, they do not require e-Signatures unless the **Include** checkbox is checked).

- For instructions to include selected actions for e-Signature requirements, see “Select the actions that require e-Signature” on page 910.

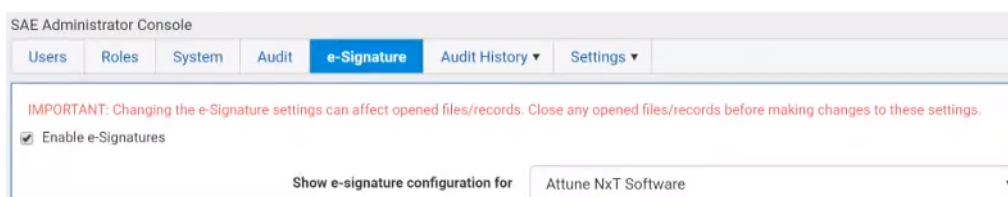
- The **e-Signature Meanings** are the stated reason for an e-Signature. There two built-in **e-Signature Meanings**:
  - **Require and Approve Experiment for Template**
  - **Review and Approve Data**

If desired, you can add custom meanings (see “Audit History tab” on page 913).

- You can set the number of signatures required for each selected action and for each required meaning in the **Number of signatures required for selected action** panel (see “Select the actions that require e-Signature” on page 910).

## Enable or disable the e-Signature function

1. In the SAE Administrator Console main screen, click the **e-Signature** tab.
2. Select **Enable e-Signature**.




---

**IMPORTANT!** Changing the e-Signature settings can affect opened files or records. Close any opened files or records before making changes to e-Signature settings.

---

3. From the **Show e-Signature configuration for** dropdown, select **Attune™ Software**.

---

**Note:** If your Attune™ Cytometric Software is configured to manage the **SAE module** for more than one application, you can configure the e-Signature settings for more than one application (for example, Attune™ Cytometric Software and SeqStudio™ Genetic Analyzer).

---

4. (Optional) Set or modify the e-Signature settings.
5. Click **Apply Settings**.

## Select the actions that require e-Signature

1. In the SAE Administrator Console main screen, click the **e-Signature** tab.
2. In the **Actions Requiring e-Signatures** panel, select each action that requires an e-Signature.

Actions Requiring Signatures

Include	Action
<input checked="" type="checkbox"/>	Export Data
<input type="checkbox"/>	Print Data
<input type="checkbox"/>	Save as Template

- For each meaning of each selected action, enter the number of e-Signatures that are required from each SAE role before that action can be performed.

Number of signatures required for selected action

Meanings	Administrator	Advanced User	User
Review and Approve Experiment for Template	0	0	0
Review and Approve Data	0	0	0

- Click **Apply Settings**.

---

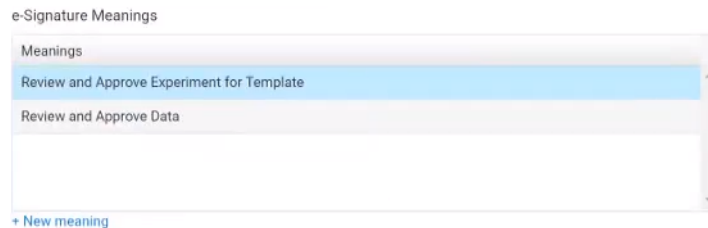
**Note:** Actions that require signatures to be completed cannot be performed unless the required signatures are provided.

---

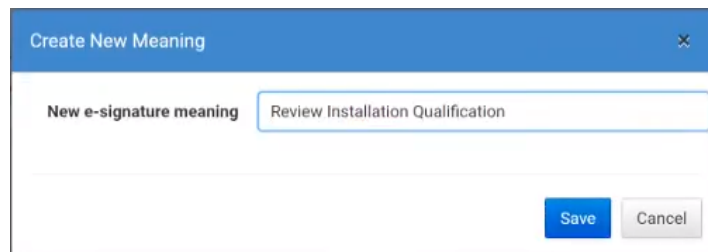
## Configure the meanings of e-Signatures

### Add an e-Signature meaning

- In the SAE Administrator Console main screen, click the **e-Signature** tab.
- In the **e-Signature Meanings** panel, click **New meaning**.



- In the **Create New Meaning** dialog, enter an e-Signature meaning in the **Name** field, then click **Save**.



- In the **e-Signature Meanings** panel, select a meaning from the **Meanings** list.
- In the **Data signed for the selected meaning** list, select the item with which to associate the meaning.

Data signed for selected meaning

Sign	Data
<input checked="" type="checkbox"/>	Tube or Plate Experiment



6. Set the actions that require e-Signatures and the number of e-Signatures that are required for that action (see “Select the actions that require e-Signature” on page 910).
7. Click **Apply Settings**.

### Delete an e-Signature meaning

1. In the SAE Administrator Console main screen, click the **e-Signature** tab.
2. In the **e-Signature Meanings** pane, select a meaning from the **Meanings** list, then click **Delete**.

---

**Note:** Default meanings cannot be deleted.

---

3. Confirm the deletion of the meaning, then click **OK**.
4. Click **Apply Settings**.

## Audit History tab

When SAE mode is active and application auditing is enabled, any change to an Experiment results in an audit transaction that is captured in the SAE console where the old and new values of each Experiment attribute is recorded.

**Audit History** tab allows users with permission to view Audit History to access the audit logs for the auditable transactions. The permission to view the entire Audit History or specific audit logs is granted in the **Roles** tab (“Roles tab” on page 886).



- Use the **Audit History dropdown** to select the audit history to view. Audit logs available for viewing are:
  - **Action Records** (“Action Records audit log” on page 914)
  - **System Configuration Records** (“System Configuration audit log” on page 917)
  - **Application Object Records** (“Application Objects Records audit log” on page 919)
  - **Instrument Run Records** (this feature is not used by the Attune™ Cytometric Software) (“Application Objects Records audit log” on page 919)

The following example shows the **Audit History** for **Action Records**.

SAE Administrator Console English ▾ | Default Administrator

Users Roles System Audit e-Signature **Audit History** Settings ▾

Enable Action Records Filtering

Date	Account Type	User Name	Full Name	Host ID	Instrument Name	Application	Action	Comment
04-Nov-2019 10:18:01 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT	Attune NxT Software	Activate Sample	A5784B83-AD55-408D-ADA2-1B69105954BC
04-Nov-2019 10:18:01 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT	Attune NxT Software	Activate Experiment	AEA9F4C-061F-4806-86B2-77A0BF2B7ECD
04-Nov-2019 10:16:01 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT	Attune NxT Software	Sign In Success	NA
04-Nov-2019 10:14:15 PST	Local	admin	admin	USEUG-6DBK5S2	Attune NxT	Attune NxT Software	Sign Out	This is a non SAE Attune user account.
04-Nov-2019 10:06:11 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT	Attune NxT Software	Enable Security	NA
04-Nov-2019 09:44:50 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	NA	SAE Administrator Console	Sign In Success	NA
04-Nov-2019 09:42:51 PST	Local	admin	admin	USEUG-6DBK5S2	Attune NxT	Attune NxT Software	Sign In Success	This is a non SAE Attune user account.

1 - 7 of 7 items

Export Report

- For instructions about how to view specific audit logs and generate printable audit records, see “View audit logs (Audit History)” on page 924.

- If desired, you can archive the audit records or set up auto archive settings in the **Settings** tab of the **SAE Administrator Console** (“Settings tab” on page 927).

Archiving audit records removes the records from the **SAE Administrator Console** and saves them in an internally specified location on the same computer on which the **SAE Administrator Console** is installed.

Archived audit records remain available for viewing from the **Settings** tab of the **SAE Administrator Console**.

For instructions about how to archive and restore audit records, see “Settings tab” on page 927.

## Action Records audit log

**Action Records** audit history shows a record of auditable actions taken by any signed in user, including signing in and out of the application, audit events on an application (such activate, duplicate, delete), running instrument maintenance functions (Startup, Shutdown, Performance Test, etc.), including custom maintenance functions, and actions taken on the SAE Administrator Console.

SAE Administrator Console English ▾ Default Administrator

Users Roles System Audit e-Signature **Audit History ▾** Settings ▾ 👤 ▾

■ Enable Action Records Filtering

Date	Account Type	User Name	Full Name	Host ID	Instrument Name	Application	Action	Comment
04-Nov-2019 10:18:01 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT	Attune NxT Software	Activate Sample	A5784B83-AD55-408D-ADA2-1B69105954BC
04-Nov-2019 10:18:01 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT	Attune NxT Software	Activate Experiment	AEAE9F4C-061F-4806-86B2-77A0BF2B7ECD
04-Nov-2019 10:16:01 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT	Attune NxT Software	Sign In Success	NA
04-Nov-2019 10:14:15 PST	Local	admin	admin	USEUG-6DBK5S2	Attune NxT	Attune NxT Software	Sign Out	This is a non SAE Attune user account.
04-Nov-2019 10:06:11 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT	Attune NxT Software	Enable Security	NA
04-Nov-2019 09:44:50 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	NA	SAE Administrator Console	Sign In Success	NA
04-Nov-2019 09:42:51 PST	Local	admin	admin	USEUG-6DBK5S2	Attune NxT	Attune NxT Software	Sign In Success	This is a non SAE Attune user account.

1 50 items per page 1 - 7 of 7 items

[Export](#) [Report](#)

- **Action Records** are always audited and cannot be turned off.
- All items in the **Action Records** are audited silently (see “Audit Mode”, “Audit Settings panel” on page 905).
- **Actions Records** contains the following columns:
  - Date
  - Account Type (Local or SAE)
  - Username
  - Full Name
  - Host ID
  - Instrument Name
  - Application Name



- Action
- Comment
- The **Host ID** is the PC name or instrument serial number.
- **Application Name** is the application for which the audit log is created.

---

**Note:** If the **SAE Administrator Console** is configured to manage the SAE module for more than one application, the **Action Records** can include audit records for the configured applications (for example, Attune™ Cytometric Software and SeqStudio™ Genetic Analyzer).

---

- If the action was taken by a non-SAE account, the **Comment** column displays “This is a non-SAE Attune™ user account.”
- If the audited action affects an application object (such as Experiment, Sample, etc.), the **Comment** column displays the GUID of the application object.

The GUID is the unique ID assigned to each application object. It is an immutable property and cannot be changed.

In the following example, the **Comments** column displays the GUID of the Sample and the Experiment listed as “Activated” in the **Action** column.

Action	Comment
Activate Sample	A5784B83-AD55-408D-ADA2-1B69105954BC
Activate Experiment	AEAE9F4C-061F-4806-86B2-77A0BF2B7ECD
Sign In Success	NA
Sign Out	This is a non SAE Attune user account.

- In addition to instrument maintenance actions that are automatically added to the **Action Records** audit history, the **Action** and **Comments** columns also include the custom maintenance functions that are added using the **Maintenance** tab of the **System Log dialog** (see “Maintenance tab” on page 800).

Action	Comment
System Maintenance Perfo...	System maintenance performed: I am a custom maintenance action too
Fluidics Filters Replaced	Fluidics Filter Replaced
Sample Syringe Replaced	Sample Syringe Replaced
Run SIP Sanitize	Instrument
Run Rinse	NA

- To filter the **Actions Records**, select the **Enable Action Records Filtering** option, then select the desired **Application** and **Action** from the corresponding dropdown list.



If desired, select or enter other filter criteria using the available controls.

Click **Search** to apply the filter criteria to the **Application Object Records**.



- Click **Report** to generate a PDF file of the log.
- Click **Export** to generate a TXT file of the log.

### Auditable actions in the SAE console

Auditable actions in the **SAE Administrator Console** are:

- Enable or disable security, audit, and e-Signature
- Sign in or out of the SAE Administrator Console
- Import or export an SAE configuration
- Install an application profile
- Archive, purge, or restore audit records
- Manual Sync with LDAP Directory



## System Configuration audit log

**System Configuration** audit history displays a record of the system security, audit, and e-Signature configuration records.

SAE Administrator Console

English | Default Administrator

Users Roles System Audit e-Signature **Audit History** Settings

Enable System Configuration Records Filtering

Changed Date	Account Ty...	Change By	Full Name	Record Type	Record Name	Action	Changed Field	Old Value	New Value
04-Nov-2019 09:59:56 PST	Local	Administrator	Default Administrator	Account	TestUser	Add	Assigned Role	NA	TestRole
				Account	TestUser	Add	First Name	NA	Test
							Last Name	NA	User
							User Name	NA	TestUser
							Password	NA	*****
							Account Type	NA	LOCAL
							Change Password at Sign In	NA	false
							Account Status	NA	ACTIVE
							Last Password Changed Date	NA	04-Nov-2019 09:59:56 PST
				04-Nov-2019 09:59:02 PST	Local	Administrator	Default Administrator	Role	TestRole
Role	TestRole	Add	Allowed Functions					NA	Export experiments (Attune NxT ...
Role	TestRole	Add	Allowed Functions					NA	Run recover sample (Attune NxT ...
Role	TestRole	Add	Allowed Functions					NA	Modify plot statistics (Attune NxT ...
Role	TestRole	Add	Allowed Functions					NA	Modify display options (Attune Nx...
Role	TestRole	Add	Allowed Functions					NA	Format plot text (Attune NxT Sof...
Role	TestRole	Add	Allowed Functions					NA	Edit text box text (Attune NxT Sof...
Role	TestRole	Add	Allowed Functions					NA	Run auto sampler calibration (Att...
Role	TestRole	Add	Allowed Functions					NA	Align objects (Attune NxT Softwa...
Role	TestRole	Add	Allowed Functions					NA	Reset baseline (Attune NxT Softw...
Role	TestRole	Add	Allowed Functions					NA	Perform overlay calculations (Att...
Role	TestRole	Add	Allowed Functions					NA	Change plot scale (Attune NxT So...
Role	TestRole	Add	Allowed Functions					NA	Change overlay plot color (Attune ...
Role	TestRole	Add	Allowed Functions					NA	Change gate type (Attune NxT So...
Role	TestRole	Add	Allowed Functions					NA	Export sample list (Attune NxT So...
Role	TestRole	Add	Allowed Functions					NA	Format statistics (Attune NxT Sof...
Role	TestRole	Add	Allowed Functions					NA	Format text box (Attune NxT Sof...
Role	TestRole	Add	Allowed Functions					NA	View system logs (Attune NxT So...
Role	TestRole	Add	Allowed Functions					NA	Import compensation (Attune Nx...
Role	TestRole	Add	Allowed Functions					NA	Run deep clean and shutdown (AT...
Role	TestRole	Add	Allowed Functions	NA	Open experiments (Attune NxT S...				

1 - 3 of 3 items

Export Report

- **System Configuration** is always audited and cannot be turned off.
- **System Configuration** audit log contains the following columns:
  - Changed Date
  - Account Type (Local or SAE)
  - Change By
  - Full Name
  - Record Type
  - Record Name
  - Action
  - Changed Field
  - Old Value
  - New Value
- The **Action** field shows the auditable action taken by the user listed in the **Change By** and **Full Name** columns.
- The **Record Type**, **Record Name**, **Changed Field**, **Old Value**, and **New Value** columns show the changes made to the system configuration, and the old and new values for the changed configuration.

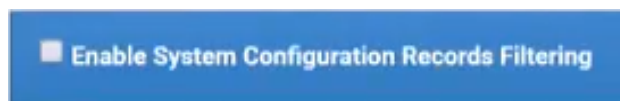
- When a change is made to the system configuration, such as add or delete users, create or assign roles, change system security settings, change audit and e-Signature configurations, it results in an audit change record.

Only the attributes that are modified are logged and the old value vs. new value is recorded in the SAE console.

In the following example, the **System Configuration** audit log shows that a new SAE Account with the account name TestUser is added to the application and the assigned role of the TestUser is TestRole.

Record Type	Record Name	Action	Changed Field	Old Value	New Value
Account	TestUser	Add	Assigned Role	NA	TestRole
Account	TestUser	Add	First Name	NA	Test
			Last Name	NA	User
			User Name	NA	TestUser
			Password	NA	*****
			Account Type	NA	LOCAL
			Change Password at Sign In	NA	false
			Account Status	NA	ACTIVE
			Last Password Changed Date	NA	04-Nov-2019 09:59:56 PST

- To filter the **System Configuration Records** audit log, select the **Enable System Configuration Records Filtering** option, then select the desired **Record type** from the corresponding dropdown list.



If desired, select or enter other filter criteria using the available controls:

- Click **Search** to apply the filter criteria to the **Application Object Records**.
- Click **Report** to generate a PDF file of the **System Configuration** audit log.
- Click **Export** to generate a TXT file of the **System Configuration** audit log.

## Application Objects Records audit log

**Application Object Records** audit history enables you to view a log of the application object records.

SAE Administrator Console English ▾ | Default Administrator

Users | Roles | System | Audit | e-Signature | **Audit History ▾** | Settings ▾

---

**Enable Application Objects Filtering**

Application	Host ID	Instrument Name	Object Type	Object Name	Last Modified Date	Last Modified By	Last Modification Reason	Last Modification Comment
Attune NxT Software	USEUG-6DBK5S2	Attune NxT	Experiment	C3E749C7-6DF1-407D-BE5D-66A...	04-Nov-2019 10:58:33 PST	TestUser	User wants to create	User wants to create
Attune NxT Software	USEUG-6DBK5S2	Attune NxT	Experiment	961A9AF3-8D4A-4053-B48A-3DE...	04-Nov-2019 10:53:36 PST	TestUser	User wants to update	User wants to update
Attune NxT Software	USEUG-6DBK5S2	Attune NxT	Experiment	BCC9253B-A92C-4CBE-8C66-DF4...	04-Nov-2019 10:49:27 PST	TestUser	User wants to update	User wants to update
Attune NxT Software	USEUG-6DBK5S2	Attune NxT	Experiment	AEA9F4C-061F-4806-86B2-77A...	04-Nov-2019 10:35:39 PST	Administrator	Calculation error.	User wants to update

1 - 4 of 4 items

---

**Data Audits (29)** | **e-Signature Records**

Changed Date	Change By	Full Name	Record Type	Record Name	Action	Changed Field	Old Value	New Value
04-Nov-2019 10:58:33 PST	TestUser	Test User	Experiment	C3E749C7-6DF1-407D-BE5...	Add	Name Notes Owner ID Owner Name Color Is Template Instrument Model Compensation Enabled Compensation Source Source Experiment ID Source Host ID	NA NA NA NA NA NA NA NA NA NA NA	New Name(1) NA F4284D40-1C03-46DA-9ADE-FC049... TestUser RGB(147, 149, 152) FALSE 4486521 FALSE NA 961A9AF3-8D4A-4053-B48A-3DEA4... USEUG-6DBK5S2
04-Nov-2019 10:58:33 PST	TestUser	Test User	Group	241C038A-B079-44FF-A73...	Add	Name Notes Owner ID Owner Name Color	NA NA NA NA NA	Group NA C3E749C7-6DF1-407D-BE5D-66A85... New Name(1) RGB(59, 114, 147)

1 - 29 of 29 items

[Report](#)

- The **Audit History** ▶ **Application Objects** tab is divided into two panels:
  - **Application objects list** (upper panel; “Application Objects list” on page 920)
  - **Data Audits** and **e-Signature Records** tabs (bottom panel; “Data Audits tab” on page 922 and “e-Signature Records tab” on page 923, respectively)

## Application Objects list

The **Application Objects** panel is the upper panel in the **Audit History** ▶ **Application Objects** tab, and it displays a list of the audited application objects.

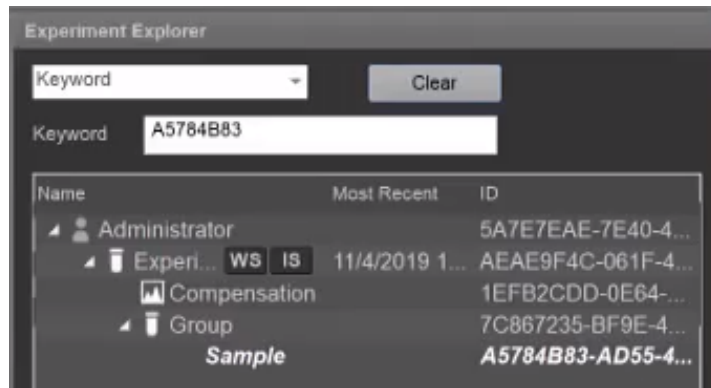
Enable Application Objects Filtering								
Application	Host ID	Instrument Name	Object Type	Object Name	Last Modified Date	Last Modified By	Last Modification Reason	Last Modification Comment
Attune NxT Software	USEUG-6DBK5S2	Attune NxT	Experiment	C3E749C7-6DF1-407D-BE5D-66A...	04-Nov-2019 10:58:33 PST	TestUser	User wants to create	User wants to create
Attune NxT Software	USEUG-6DBK5S2	Attune NxT	Experiment	961A9AF3-8D4A-4053-B48A-3DE...	04-Nov-2019 10:53:36 PST	TestUser	User wants to update	User wants to update
Attune NxT Software	USEUG-6DBK5S2	Attune NxT	Experiment	BCC9253B-A92C-4CBE-8C66-DF4...	04-Nov-2019 10:49:27 PST	TestUser	User wants to update	User wants to update
Attune NxT Software	USEUG-6DBK5S2	Attune NxT	Experiment	AEA9F4C-061F-4806-86B2-77A...	04-Nov-2019 10:35:39 PST	Administrator	Calculation error.	User wants to update

- Application objects are auditable items such as plate setups, templates, or other items that you create in an application.
- Examples of application object types in the Attune™ Cytometric Software are **Plate Experiment**, **Tube Experiment**, **Template**, **Filter Configuration**, and **Filter Definition**.
- When an application object is created, updated, or deleted, an audit record is generated on the SAE console. An application object can have multiple data audit records and multiple e-Signature records.
- The **Application Objects** list has the following columns:
  - Application
  - Host ID
  - Instrument Name
  - Object Type
  - Object Name
  - Last Modified Date
  - Last Modified By
  - Last Modification Reason
  - Last Modification Comment
- The **Host ID** is the PC name or instrument serial number.
- The **Object Name** is the GUID of the object. The GUID is a unique ID assigned each an application object. It is an immutable property and cannot be changed.
- Each application object record has a corresponding audit log consisting of all subapplication object components with their attribute value change record (old value vs. new value).  
If you have permission to view **Application Object Records** in **Audit History**, you can view the audit records in the **Data Audits** tab (“Data Audits tab” on page 922).
- If the **e-Signature** feature is enabled in the **e-Signature** tab (“e-Signature tab” on page 909) and you have permission to view **Application Object Records** in **Audit History**, you can view the e-Signature log for the selected application object in the **e-Signature Records** tab (“e-Signature Records tab” on page 923).
- If an application object is deleted, the deleted object appears with a strike-through in the SAE server.

In the following example, the second experiment from the top is deleted. As a result, the GUID of the deleted experiment is shown with a line over it (strike-through).

Application	Host ID	Instrument Name	Object Type	Object Name	Last Modified
Attune NxT Software	USEUG-6DBK5S2	Attune NxT	Experiment	AE48102E-79FF-4A93-ABF6-8508...	04-Nov-
Attune NxT Software	USEUG-6DBK5S2	Attune NxT	Experiment	961A9AF3-804A-4053-B48A-3DE...	04-Nov-
Attune NxT Software	USEUG-6DBK5S2	Attune NxT	Experiment	C3E749C7-6DF1-407D-8E5D-66A...	04-Nov-

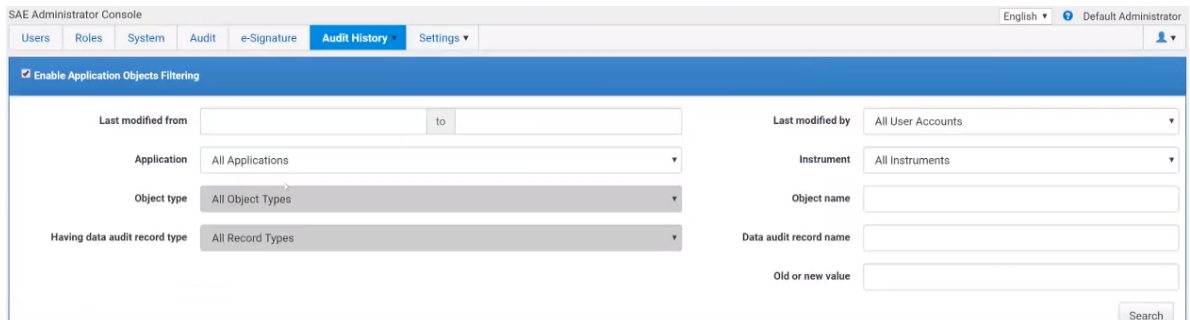
**Note:** When the Attune™ Cytometric Software is in the SAE mode, you can view the GUID of application objects in the **Experiment Explorer**, and perform a **Keyword** search using the objects GUID (see “Display ID” on page 299).



- To view a list of items that are audited for your application, select **Enable Application Object Records Filtering**.

In the **Filtering** panel, select Attune™ Cytometric Software for the **Application**, then select or enter the desired filter criteria using the available controls.

Click **Search** to apply the filter criteria to the **Application Object Records**.



## Data Audits tab

The **Data Audits** tab in the bottom panel contains the audit log corresponding to the selected application object.

Data Audits (29)		e-Signature Records							
Changed Date	Change By	Full Name	Record Type	Record Name	Action	Changed Field	Old Value	New Value	
04-Nov-2019 10:58:33 PST	TestUser	Test User	Experiment	C3E749C7-6DF1-407D-BE5...	Add	Name	NA	New Name(1)	
						Notes	NA	NA	
						Owner ID	NA	F4284D40-1C03-46DA-9ADE-FC049...	
						Owner Name	NA	TestUser	
						Color	NA	RGB(147, 149, 152)	
						Is Template	NA	FALSE	
						Instrument Model	NA	4486521	
						Compensation Enabled	NA	FALSE	
						Compensation Source	NA	NA	
						Source Experiment ID	NA	961A9AF3-8D4A-4053-B48A-3DEA4...	
						Source Host ID	NA	USEUG-6DBK5S2	
04-Nov-2019 10:58:33 PST	TestUser	Test User	Group	241C038A-B079-44FF-A73...	Add	Name	NA	Group	
						Notes	NA	NA	
						Owner ID	NA	C3E749C7-6DF1-407D-BE5D-66A85...	
						Owner Name	NA	New Name(1)	
						Color	NA	RGB(59, 114, 147)	
04-Nov-2019 10:58:33 PST	TestUser	Test User	Sample	AC99BB63-B6ED-400A-AED...	Add	Name	NA	Sample(1)	

50 Items per page 1 - 29 of 29 items

- The **Data Audits** tab has the following columns:
  - Changed Date
  - Change By
  - Full Name
  - Record Type
  - Record Name
  - Action
  - Changed Field
  - Old Value
  - New Value





- The **Changed Field**, **Old Value**, and **New Value** columns show the subapplication object components of an application object, and the new and old values of its attributes. When an application object is first created, each of these attributes displays **NA** for the **Old Value**.

When any of these attributes are changed after the object is created, it results in an audit change record. Only the attributes that are modified are logged and the old value versus the new value are recorded in the SAE console.

The following example shows the **Changed Field**, **Old Value**, and **New Value** columns for a **Plate Record**.

Changed Field	Old Value	New Value
Name	NA	Plate(2)
Notes	NA	NA
Owner ID	NA	B817D128-7609-42BD-9831-07412...
Owner Name	NA	Administrator
Description	NA	NA
Barcode (Plate ID)	NA	NA
Is Template	NA	FALSE
Instrument Model	NA	4486521
Run Orientation	NA	Horizontal
Collection Mode	NA	Only Wells - 1,2,3,4,5,6,7,13,14,15,1...
Collection Wells	NA	A1, A10, A11, A12, A2, A3, A4, A5, ...
Source Plate ID	NA	C2DEF134-4D36-4F0B-B771-A2C9...
Source Host ID	NA	USEUG-6DBK5S2

- Click **Report** to generate a PDF file of the **Data Audits** log. **Export** is not supported for the **Data Audits** log.

### e-Signature Records tab

The **e-Signature Records** tab in the bottom panel displays the record of e-Signatures and the status of the signatures for the selected application object.

Data Audits (55)		e-Signature Records (1)			
Signature Date	Signature Meaning	Signed By	Full Name	Role	Status
04-Nov-2019 11:20:27 PST	Review and Approve Data	Administrator	Default Administrator	Administrator	CURRENT

10 items per page | 1 - 1 of 1 items

- The **e-Signature Records** tab has the following columns:
  - Signature Date
  - Signature Meaning
  - Signed By
  - Full Name
  - Role
  - Status (Current or Obsolete)

- Signatures represent the state of the Experiment at the time of signing. If any modifications are made to the Experiment after signing, the existing signature becomes obsolete.

Data Audits (55)		e-Signature Records (1)			
Signature Date	Signature Meaning	Signed By	Full Name	Role	Status
04-Nov-2019 11:20:27 PST	Review and Approve Data	Administrator	Default Administrator	Administrator	OBSOLETE

- When the signatures become obsolete, actions that require signatures cannot be performed unless new signatures are provided.
- Click **Report** to generate a PDF file of the **e-Signature Records** log.

## Instrument Run Records audit log

**Instrument Run Records** audit history shows a record of the instrument run, including a summary of the run, the application objects used and the actions performed in the run, any changes made during the run, and a list outputs at the completion of the run.

---

**IMPORTANT!** **Instrument Run Records** audit log is not used by the Attune™ Cytometric Software. Therefore, the **Instrument Run Records** audit history is not available for the Attune™ Cytometric Software application.

---

## View audit logs (Audit History)

To view audit logs in the **Audit History** tab, you need to have permission to view **Audit History** or specific audit logs, which is granted in the **Roles** tab (“Roles tab” on page 886).



## View the Action Records audit log

All items in the **Action Records** log are audited silently.

1. In the **SAE Administrator Console** main screen, click the **Audit History** tab.
2. Select **Action Records** to view a log of the specified audit events.
3. To view a list of items that are audited for the application:
  - a. Select **Enable Action Records Filtering**.
  - b. Select the application from the **Application** list.
  - c. In the **Action** field, click the dropdown arrow to view the list of auditable actions.



4. (Optional) Perform the following actions:
  - Specify other filtering settings.
  - Click **Report** to generate a PDF file of the log.
  - Click **Export** to generate a TXT file of the log.

### View the System Configuration audit log

1. In the **SAE Administrator Console** main screen, click the **Audit History** tab.
2. **Select System Configuration** to view a log of the System Security, Audit, and e-Signature configuration records.
3. To view a list of items that are audited:
  - a. Select **Enable System Configuration Records Filtering**.
  - b. In the **Record Type** field, click the dropdown arrow to view the list of auditable system configuration objects.
4. (Optional) Perform the following actions:
  - Specify other filtering settings.
  - Click **Report** to generate a PDF file of the log.
  - Click **Export** to generate a TXT file of the log.

### View the Application Object Records audit log

Application objects are auditable items such as plate setups, templates, or other items that you create in an application.

1. In the **SAE Administrator Console** main screen, click the **Audit History** tab.
2. Select **Application Object Records**.
3. To view a list of items that are audited for the application:
  - a. Select **Enable Application Object Records Filtering**.
  - b. Select the application from the **Application** list.
  - c. In the **Having data audit record type** field, click the dropdown arrow to view the list of auditable objects.
4. (Optional) Perform the following actions:
  - Specify other filtering settings.
  - Click **Report** to generate a PDF file of the log.
  - **Export** is not supported for the **Application Object Records** audit log.



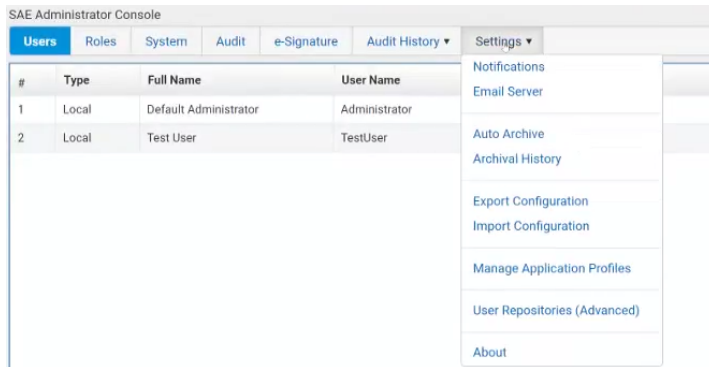
## View the Instrument Run Records audit log

Audit records are listed when a run is complete.

1. In the **SAE Administrator Console** main screen, click the **Audit History** tab.
2. Select **Instrument Run Records**, then select the tab for the specific audit log you want to view:
  - Run Summary
  - Application Objects
  - Action Records
  - Data Audit Records
  - Run Completion Outputs
3. To limit the records that are displayed, select **Enable Instrument Run Records Filtering**.
4. *(Optional)* Perform the following actions:
  - Specify other filtering settings.
  - Click **Report** to generate a PDF file of the log.
  - **Export** is not supported for the **Instrument Run Records** audit log.

## Settings tab

**Settings** tab allows users that have permission to modify **SAE Administrator Console** to configure optional and advanced SAE Console settings, which are available as a dropdown list from the **Settings** tab.



Using the **Settings** dropdown, you can view or configure the following options:

- **Notifications:** Enables you to specify when and how to be notified when specified events occur in the SAE (see “Set up SAE messaging notifications” on page 928).
- **Email Server:** Enables you to configure the SMTP server so that the SAE Administrator Console can send email notifications (see “Configure the SMTP server for email notifications” on page 929).
- **Auto Archive:** Enables you to archive audit records automatically (see “Auto archive audit records” on page 929).
- **Archival History:** Enables you to:
  - Archive audit records manually (see “Archive audit records manually” on page 930).
  - View or export archived audit records (see “View or export archived audit records” on page 930).
  - Restore archived audit records (see “Restore archived audit records” on page 931).
- **Export Configuration:** Enables you to export user, system security, audit, and e-Signature settings from the **SAE Administrator Console** (see “Export configuration” on page 932).
- **Import Configuration:** Enables you to import settings from another installation of the **SAE Administrator Console** (see “Import configuration” on page 932).
- **Manage Application Profiles:** Enables you to install a profile for the application for which you want to configure the SAE module (see “Set up the SAE administrator console with application profiles” on page 933).
- **User Repositories (Advanced):** Enables you to configure user repositories for SAE or external account access (see “Configure user repositories” on page 933).
- **About:** Displays the **SAE Administrator Console** software version.

## Set up SAE messaging notifications

You can specify when and how to be notified when specified events occur in the SAE. You have the option to be notified at sign in, to be notified by email when the event of interest occurs, or both.

---

**IMPORTANT!** You **must** configure the SMTP server before you set up email notifications (“Configure the SMTP server for email notifications” on page 929).

---

1. In the SAE Administrator Console main screen, click the **Settings** tab, then select **Notifications**.

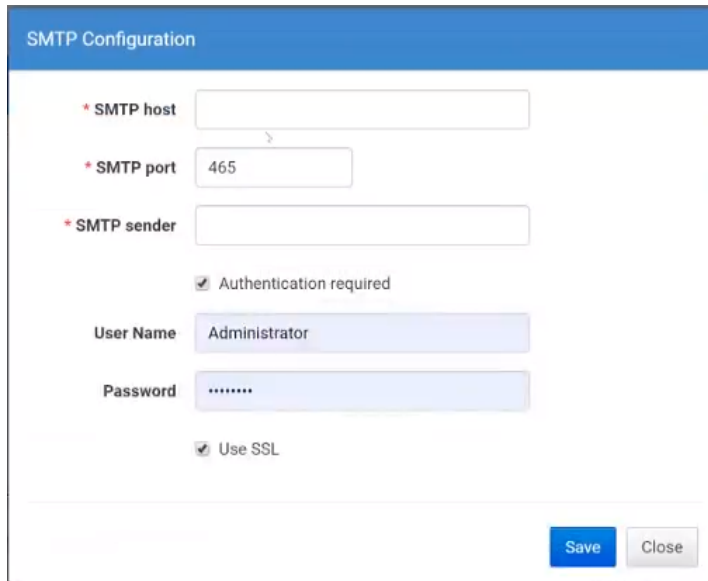
Event	Notify at Administrator Sign In	Notify by Email	Email Address
Security enabled	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Separate multiple emails (max. 5) using comma
Security disabled	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Separate multiple emails (max. 5) using comma
User did not enter correct password	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	Separate multiple emails (max. 5) using comma
User account suspended	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	Separate multiple emails (max. 5) using comma
User session timeout	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Separate multiple emails (max. 5) using comma
Role deleted	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Separate multiple emails (max. 5) using comma

2. In the **Edit Notification Settings** dialog, select **Notify at Administrator sign in** for the events of interest.
3. (Optional) Select **Notify by Email**, then specify an email address.
4. Click **Save**.
  - If you have selected **Notify at Administrator sign in**, you will be notified at each sign in if an event of interest occurs while you were signed out.
  - If you have selected **Notify by Email**, an email will be sent to you when an event of interest occurs.

## Configure the SMTP server for email notifications

Before you set up email notifications (“Set up SAE messaging notifications” on page 928), you must configure the SMTP server so that the SAE Administrator Console can send email notifications.

1. In the SAE Administrator Console main screen, click the **Settings** tab, then select **Email Server**.



The screenshot shows the 'SMTP Configuration' dialog box. It has a blue header with the text 'SMTP Configuration'. Below the header are several input fields and checkboxes. The fields are: '\* SMTP host' (empty), '\* SMTP port' (465), '\* SMTP sender' (empty), 'User Name' (Administrator), and 'Password' (masked with dots). There are two checkboxes: 'Authentication required' (checked) and 'Use SSL' (checked). At the bottom right of the dialog are two buttons: 'Save' (blue) and 'Close' (grey).

2. In the **SMTP Configuration** dialog, enter the following:

- **SMTP host, SMTP port, and SMTP sender**

---

**Note:** Select **Authentication required** if the SMTP server requires authentication.

---

- **Username and Password**

---

**Note:** Select **Use SSL** if the SMTP server requires an encrypted channel connection.

---

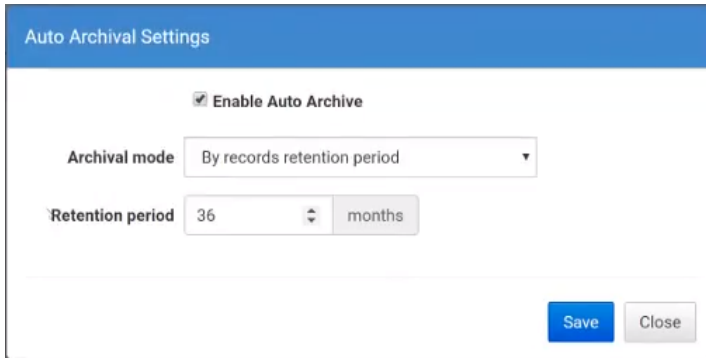
3. Click **Save**.

## Auto archive audit records

Archiving audit records removes the records from the SAE Administrator Console and saves them in an internally specified location on the same computer on which the SAE Administrator Console is installed.

1. In the SAE Administrator Console main screen, click the **Settings** tab.
2. Select **Auto Archive**.

3. In the **Auto Archival Settings** dialog, select **Enable Auto Archive**.



4. Select the **Archival mode** and associated settings:

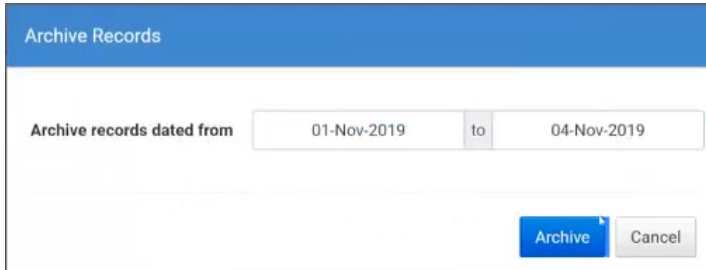
- By number of records or retention period
- By number of records
- By retention period

5. Click **Save**.

The software periodically checks the audit record status and archives when the specified archive conditions are met.

## Archive audit records manually

1. In the SAE Administrator Console main screen, click the **Settings** tab.
2. Select **Archival History**.
3. Click **Ad-hoc Archive**, select the start and end dates, then click **Archive**.



---

**Note:** Archived audit records (automatic or manual) are accessible for viewing in the SAE Administrator Console (see “View or export archived audit records” on page 930).

---

## View or export archived audit records

1. In the SAE Administrator Console main screen, click the **Settings** tab.
2. Select **Archival History**.

The list of archived records is displayed in the **Settings tab** ▶ **Archival history** screen.





Run Date	Run Duration	Records Dated From	Records Dated To	Records Count	Server File Path
04-Nov-2019 11:35:37 PST	307 ms	01-Nov-2019 00:00:00 PDT	04-Nov-2019 00:00:00 PST	Action records : 49 System configuration records : 13 Application objects records : 5 Instrument run records : 0	C:\Program Files\ThermoFisher\SAEAdmin\automated_archivals\04-Nov-2019 11:35:37-571 PST

- To view a record, select the record from the list, then click **View Archived Records**. Audit log for the selected record is displayed, arranged in **Action Records, System Configuration, Application Object Records, and Instrument Run Records** tabs.

Date	Account	User Name	Full Name	Host ID	Application	Action	Comment
04-Nov-2019 09:42:51 PST	Local	admin	admin	USEUG-6DBK5S2	Attune NxT Software	Sign In Success	This is a non SAE Attune user account.
04-Nov-2019 09:44:50 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	SAE Administrator Co...	Sign In Success	NA
04-Nov-2019 10:06:11 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT Software	Enable Security	NA
04-Nov-2019 10:14:15 PST	Local	admin	admin	USEUG-6DBK5S2	Attune NxT Software	Sign Out	This is a non SAE Attune user account.
04-Nov-2019 10:16:01 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT Software	Sign In Success	NA
04-Nov-2019 10:18:01 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT Software	Activate Experiment	AEAE9F4C-061F-4806-86B2-77A0BF2B7ECD
04-Nov-2019 10:18:01 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT Software	Activate Sample	A5784B83-AD55-408D-ADA2-1B69105954BC
04-Nov-2019 10:25:19 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT Software	Activate Experiment	AEAE9F4C-061F-4806-86B2-77A0BF2B7ECD
04-Nov-2019 10:25:19 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT Software	Activate Sample	A5784B83-AD55-408D-ADA2-1B69105954BC
04-Nov-2019 10:38:02 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT Software	Sign In Success	Application unlocked
04-Nov-2019 10:38:02 PST	Local	Administrator	Administrator	USEUG-6DBK5S2	Attune NxT Software	Sign In Success	Application unlocked
04-Nov-2019 10:38:35 PST	Local	Administrator	Default Administrator	USEUG-6DBK5S2	Attune NxT Software	Sign Out	NA
04-Nov-2019 10:38:47 PST	Local	TestUser	Test User	USEUG-6DBK5S2	Attune NxT Software	Sign In Success	NA
04-Nov-2019 10:40:42 PST	Local	TestUser	Test User	USEUG-6DBK5S2	Attune NxT Software	Activate Sample	A776C4FA-6358-40D9-96CD-ED1E3300CAE0
04-Nov-2019 10:40:42 PST	Local	TestUser	Test User	USEUG-6DBK5S2	Attune NxT Software	Activate Experiment	BCC9253B-A92C-4CBE-8C66-DF4761601EB7
04-Nov-2019 10:43:33 PST	Local	TestUser	Test User	USEUG-6DBK5S2	Attune NxT Software	Sign Out	NA
04-Nov-2019 10:44:52 PST	Local	TestUser	Test User	USEUG-6DBK5S2	Attune NxT Software	Sign In Success	NA
04-Nov-2019 10:45:23 PST	Local	TestUser	Test User	USEUG-6DBK5S2	Attune NxT Software	Activate Experiment	File Tampering Detected: A776C4FA-6358-40D9-9
04-Nov-2019 10:46:25 PST	Local	TestUser	Test User	USEUG-6DBK5S2	Attune NxT Software	Activate Experiment	File Tampering Detected: A776C4FA-6358-40D9-9
04-Nov-2019 10:46:42 PST	Local	TestUser	Test User	USEUG-6DBK5S2	Attune NxT Software	Activate Experiment	File Tampering Detected: A776C4FA-6358-40D9-9

1 2 3 20 items per page

[Back to Archival History](#)

- Click **Back to Archival History** to view the list of archived records again.
- To export a record, select the record from the **Archival history** list, then click **Export**.

## Restore archived audit records

- In the **SAE Administrator Console** main screen, click the **Settings** tab.
- Select **Archival History**.

The list of archived records is displayed in the **Settings tab** ▶ **Archival history** screen.

SAE Administrator Console						
Users	Roles	System	Audit	e-Signature	Audit History	Settings
Run Date	Run Duration	Records Dated From	Records Dated To	Records Count	Server File Path	
04-Nov-2019 11:35:37 PST	307 ms	01-Nov-2019 00:00:00 PDT	04-Nov-2019 00:00:00 PST	Action records : 49 System configuration records : 13 Application objects records : 5 Instrument run records : 0	C:\Program Files\ThermoFisher\SAEAdmin\automated_archivals\04-Nov-2019 11-35-37-571 PST	

- To restore a record, select the record from the list, then click **Restore**.
- To restore a ZIP file that was exported from the archival history, click **Restore (upload)**, then select the ZIP file.

## Export configuration

- In the SAE Administrator Console main screen, click the **Settings** tab.
- Select **Export Configuration**.
- In the **Export Configuration** dialog, select an export option:

Setting	Exports
<b>All</b>	SAE settings and SAE user accounts
<b>Custom Users &amp; Roles</b>	<ul style="list-style-type: none"> <li>SAE user accounts with Active status</li> <li>SAE roles and their associated permissions</li> </ul>
<b>Custom System &amp; Roles</b>	<ul style="list-style-type: none"> <li>SAE settings</li> <li>SAE roles and their associated permissions</li> </ul>

- Click **Export**.  
The exported file (DAT format) downloads to the default location of the computer.

## Import configuration

- In the SAE Administrator Console main screen, click the **Settings** tab.
- Select **Import Configuration**.
- Click **Choose File** to select the DAT file with the desired configuration settings.
- Select an import option.
- Click **Import**.
- If imported user accounts exist in the SAE Administrator Console, click **Skip** or **Overwrite** for each user account, then click **Confirm and Import**.



## Set up the SAE administrator console with application profiles

An application profile contains default SAE Administrator Console settings for an application (such as the Attune™ Cytometric Software, SeqStudio™ Genetic Analyzer, etc.).

Before you can use the SAE Administrator Console to configure the SAE module for an application, you must install a profile for the application.

1. In the SAE Administrator Console main screen, click the **Settings** tab.
2. Select **Manage Application Profiles**.
3. Click **Install Application Profile**, then select the DAT file for the application profile.
4. Click **Verify Data File**.
5. Select **Install new application**, then click **Install**.

The application name and settings are added to the SAE Administrator Console.

## Configure user repositories

SAE user account information is stored in a “user repository”. The SAE Administrator Console allows you to use **Internal**, **External LDAP**, or **Federated Repositories** (see “User repository overview” on page 936).

1. In the SAE Administrator Console main screen, click the **Settings** tab.
2. Select **User Repositories (Advanced™)**.
3. Select the **User repository definition**:

Setting	Exports
<b>Internal User Repository</b>	Allows SAE user accounts to sign in
<b>External LDAP User Repository</b>	Allows external user accounts to sign in
<b>Federated Repositories</b>	Allows SAE user accounts or LDAP accounts to sign in

**Note:** For more information on SAE Administrator Console user repositories, see “User repository overview” on page 936.

4. If you selected **External LDAP User Repository** or **Federated Repositories**, click **Next**, then enter the required information (see “User repository settings” on page 934).
5. Click the **Users** tab to display the list of accounts added to the SAE Administrator Console.
  - New LDAP accounts are listed as **External**, and **Role** is set to the default specified during account mapping. If no default was specified, accounts are set to **No Privileges Role**.
  - SAE user accounts that were previously created in the SAE Administrator Console are listed as **Local**.
  - If you selected LDAP, the **Status** for all accounts except for the default SAE Administrator account is set to **Inactive**.

6. Click **Test Connection** to synchronize the new accounts with the LDAP server.  
The SAE server also periodically syncs the LDAP accounts with the LDAP server if changes are made to the User repository definition or any setting on the LDAP server.
7. If needed, edit the user accounts to assign roles.

## User repository settings

The following table lists the **External LDAP User Repository** and **Federated Repositories** settings.

Setting	Description
<b>LDAP Server Configuration</b>	
Host name, Port, and Use ssl	LDAP server name or IP address, port, and interface protocol.
Bind distinguished name, Bind password, Base distinguished name	LDAP server attributes required for access.
<b>User Account Mapping</b>	
Directory type	LDAP server configuration. Click <b>Set Defaults</b> after you select the <b>Directory type</b> to display typical default parameters for mapping to an LDAP system.
Username	Parameter that maps to the username in the LDAP system.
Default role assignment	The SAE role that will be assigned to all user accounts. You can change the role after the user accounts are imported into the SAE Administrator Console.
Username and other settings	Parameters that correspond to the username and other fields in the LDAP system.
<b>Authentication verification</b>	
Username and Password	LDAP username and password.

## Sign-in with LDAP or federated user repositories

The following table lists the required credentials for User or Administrator sign in to the SAE Administrator Console with LDAP or federated user repositories.

User repository	User signs in with	Administrator signs in with
Internal	Internal (local) account: Username and password created in the SAE Administrator Console	<ul style="list-style-type: none"> <li>• Username and password for the default SAE Administrator user account</li> <li>• Any SAE user account that has been assigned the SAE role of Administrator</li> </ul>



(continued)

User repository	User signs in with	Administrator signs in with
<b>External</b>	External account: Username and password created in the LDAP user management system.  <b>Note:</b> Local accounts are set to <b>Inactive</b> .	<ul style="list-style-type: none"> <li>• Username (with local/prefix) and password for the default SAE Administrator user account Example: local/Administrator</li> <li>• Any external account that has been assigned the SAE role of Administrator</li> </ul>
<b>Federated</b>	The account type that they are assigned: <ul style="list-style-type: none"> <li>• External account</li> <li>• Internal (local) account (with local/prefix)</li> </ul> Example: local/Username	<ul style="list-style-type: none"> <li>• Username (with local/prefix) and password for the default SAE Administrator user account Example: local/Administrator</li> <li>• Any external account that has been assigned the SAE role of Administrator</li> </ul>

## User repository overview

SAE user account information is stored in a **user repository**.

The SAE Administrator Console provides the following options for user repositories:

- **Internal:** Allows only SAE user accounts to sign in to an application. SAE user accounts are referred to as **local** accounts in the SAE Administrator Console.
  - SAE user accounts are created in the SAE Administrator Console and are identified as **local** in the **Users** tab.
  - User authentication is based on the accounts that are listed in the **Users** tab and the SAE settings that are specified in the **System** tab.
  - User permissions are determined by the roles that are configured in the SAE Administrator Console.
- **External LDAP:** Enables LDAP based authentication with an LDAP directory. Allows only **external** user accounts to sign in to an application.
  - User accounts are created in an LDAP (Lightweight Directory Access Protocol) user management system and are identified as **external** in the SAE Administrator Console **Users** tab.
  - User authentication is based on the accounts that are listed in the SAE Administrator Console **Users** tab and the external LDAP user repository. The SAE settings that are specified in the SAE Administrator Console **System** tab are not used.
  - User permissions are determined by the roles that are configured in the SAE Administrator Console.
  - All local user accounts except the default **Administrator** account are set to **Inactive**.
  - Passwords cannot be changed in the SAE Administrator Console.
- **Federated:** Allows internal (**local**) and **external** account sign-in to an application.
  - User accounts are created in the SAE Administrator Console or in an LDAP user management system.
  - User authentication is based on the respective internal or LDAP user repository.



## Overton's cumulative statistics

The Overton Subtraction method determines the positive result by subtracting a control histogram from the sample histogram on a channel-by-channel basis. Any negative differences (where the control has a greater number of cells in a channel compared to the sample) are added to the positive differences in the lower channels.

The percent positive is then calculated as the sum of the positive differences for all channels divided by the total number of events.

For a complete description of the Overton Subtraction method, refer to “Modified Histogram Subtraction Technique for Analysis of Flow Cytometry Data W. Roy Overton Cytometry 9:619-626 (1988)”.

## Kolmogorov-Smirnov test

In statistics, the Kolmogorov–Smirnov test (K–S test) is a nonparametric test for the equality of continuous, one-dimensional probability distributions that can be used to compare a sample with a reference probability distribution (one-sample K–S test) or to compare two samples (two-sample K–S test).

The Kolmogorov–Smirnov statistic quantifies a distance between the empirical distribution function of the sample and the cumulative distribution function of the reference distribution, or between the empirical distribution functions of two samples. The null distribution of this statistic is calculated under the null hypothesis that the samples are drawn from the same distribution (in the two-sample case) or that the sample is drawn from the reference distribution (in the one-sample case). In each case, the distributions considered under the null hypothesis are continuous distributions but are otherwise unrestricted.

The two-sample KS test is one of the most useful and general nonparametric methods for comparing two samples, as it is sensitive to differences in both location and shape of the empirical cumulative distribution functions of the two samples.



## Hyperlog™

Compensated flow cytometry data frequently contains negative values due to compensation, and cell populations do occur which have low means and normal distributions. Logarithmic transformations cannot properly handle negative values, and poorly display normally distributed cell types. Alternative transformations addressing this issue include the log-linear hybrid transformations and Hyperlog™.

The Hyperlog™ transform is a log-like transform that admits negative, zero, and positive values. The transform is a hybrid type of transform specifically designed for compensated data. One of its parameters allows it to smoothly transition from a logarithmic to linear type of transform that is ideal for compensated data (Cytometry A., 2005, 64:34-42).

## Linear and log color binning reference

The **Linear mode** bins data by assigning a color index for each density pixel linearly such that each increment is determined by dividing the range ( $Z_{\max} - Z_{\min}$ ) by the number of color steps.

The **Log mode** bins data by assigning a color index for each density pixel logarithmically such that each increment is determined by dividing the logarithmic range ( $\log Z_{\max} - \log Z_{\min}$ ) by the number of color steps. The index is then determined as  $\text{int}((\log Z_{\text{val}} - \log Z_{\min}) / \text{increment})$ .





# Safety



**WARNING! GENERAL SAFETY.** Using this product in a manner not specified in the user documentation may result in personal injury or damage to the instrument or device. Ensure that anyone using this product has received instructions in general safety practices for laboratories and the safety information provided in this document.

- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
- Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs) and use appropriate personal protective equipment (gloves, gowns, eye protection, and so on). To obtain SDSs, visit [thermofisher.com/support](https://www.thermofisher.com/support).

## Chemical safety



**WARNING! GENERAL CHEMICAL HANDLING.** To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below. Consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see the "Documentation and Support" section in this document.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with sufficient ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer cleanup procedures as recommended in the SDS.
- Handle chemical wastes in a fume hood.
- Ensure use of primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- After emptying a waste container, seal it with the cap provided.
- Characterize (by analysis if needed) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
- **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.



**WARNING! HAZARDOUS WASTE (from instruments).** Waste produced by the instrument is potentially hazardous. Follow the guidelines noted in the preceding General Chemical Handling warning.



**WARNING! 4L Reagent and Waste Bottle Safety.** Four-liter reagent and waste bottles can crack and leak. Each 4-liter bottle should be secured in a low-density polyethylene safety container with the cover fastened and the handles locked in the upright position.

## Biological hazard safety



**WARNING! Potential Biohazard.** Depending on the samples used on this instrument, the surface may be considered a biohazard. Use appropriate decontamination methods when working with biohazards.



**WARNING! BIOHAZARD.** Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Conduct all work in properly equipped facilities with the appropriate safety equipment (for example, physical containment devices). Safety equipment can also include items for personal protection, such as gloves, coats, gowns, shoe covers, boots, respirators, face shields, safety glasses, or goggles. Individuals should be trained according to applicable regulatory and company/ institution requirements before working with potentially biohazardous materials. Follow all applicable local, state/provincial, and/or national regulations. The following references provide general guidelines when handling biological samples in laboratory environment.

- U.S. Department of Health and Human Services, *Biosafety in Microbiological and Biomedical Laboratories (BMBL)*, 6th Edition, HHS Publication No. (CDC) 300859, Revised June 2020  
[cdc.gov/labs/bmbi](https://www.cdc.gov/labs/bmbi)
- Laboratory biosafety manual, fourth edition. Geneva: World Health Organization; 2020 (Laboratory biosafety manual, fourth edition and associated monographs)  
[who.int/publications/i/item/9789240011311](https://www.who.int/publications/i/item/9789240011311)

# Documentation and support

## Customer and technical support

Visit [thermofisher.com/support](https://www.thermofisher.com/support) for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
  - Product FAQs
  - Software, patches, and updates
  - Training for many applications and instruments
- Order and web support
- Product documentation
  - User guides, manuals, and protocols
  - Certificates of Analysis
  - Safety Data Sheets (SDSs; also known as MSDSs)

---

**Note:** For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

---

## Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](https://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have questions, contact Life Technologies at [www.thermofisher.com/support](https://www.thermofisher.com/support).



