

# Applied Biosystems 3500/3500xL Genetic Analyzer User Guide

User Guide

For Research Use Only. Not for use in diagnostic procedures.



#### For Research Use Only. Not for use in diagnostic procedures.

© Copyright 2009, 2010 Life Technologies Corporation. All rights reserved.

Information in this document is subject to change without notice.

APPLIED BIOSYSTEMS DISCLAIMS ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE. TO THE FULLEST EXTENT ALLOWED BY LAW, IN NO EVENT SHALL APPLIED BIOSYSTEMS BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF, WHETHER OR NOT FORESEEABLE AND WHETHER OR NOT APPLIED BIOSYSTEMS IS ADVISED OF THE POSSIBILITY OF SUCH DAMAGES.

#### NOTICE TO PURCHASER:

The purchase price of this Instrument includes a grant of a limited, non-transferable license under U.S. patents and method claims of its foreign counterparts, and element clams of its foreign counterparts, to use this particular instrument for electrophoresis methods employing fluorescence as a means of detection. No other licenses or rights are hereby conveyed either expressly, by implication, or estoppel including, but not limited to, any claims to a composition.

This instrument incorporates technology subject to one or more patents licensed from Hitachi, Ltd. as well as patents and patented technology owned by or under control of Applied Biosystems.

This instrument is Authorized for use in DNA sequencing and fragments analysis only. This Authorization is included in the purchase price of the instrument and corresponds to the up-front fee component of a license under process claims of U.S. patents and under all process claims for DNA sequence and fragment analysis of U.S. patents now or hereafter owned or licensable by Applied Biosystems for which an Authorization is required, and under corresponding process claims in foreign counterparts of the foregoing for which an Authorization is required. The running royalty component of licenses may be purchased from Applied Biosystems or obtained by using Authorized reagents purchased from Authorized suppliers in accordance with the label rights accompanying such reagents. Purchase of this instrument does not itself convey to the purchaser a complete license or right to perform the above processes. This instrument is also licensed under U.S. patents and apparatus and system claims in foreign counterparts thereof. No rights are granted expressly, by implication, or by estoppel under composition claims or under other process or system claims owned licensable by Applied Biosystems. For more information regarding licenses, please contact the Director of Outlicensing at Applied Biosystems, 850 Lincoln Centre Drive, Foster City. California 94404, USA.

#### TRADEMARKS:

Trademarks of Life Technologies Corporation and its affiliated companies: Applied Biosystems<sup>®</sup>, AB (Design)<sup>®</sup>, AmpFtSTR<sup>®</sup>, BigDye<sup>®</sup>, BigDye Xterminator<sup>™</sup>, COfiler<sup>®</sup>, GeneMapper<sup>®</sup>, GeneScan<sup>™</sup>, Hi-Di<sup>™</sup>, Identifiler<sup>®</sup>, KB<sup>™</sup>, LIZ<sup>®</sup>, MicroSeq<sup>®</sup>, Minifiler<sup>™</sup>, POP-4<sup>™</sup>, POP-6<sup>™</sup>, POP-7<sup>™</sup>, Profiler<sup>®</sup>, Profiler Plus<sup>®</sup>, SEfiler Plus<sup>™</sup>, SeqScape<sup>®</sup>, SGM Plus<sup>®</sup>, SNaPshot<sup>®</sup>, Variant Reporter<sup>™</sup>, Yfiler<sup>®</sup>

AFLP is a registered trademark of Keygene N.V. MLPA is a registered trademark of MRC-Holland.

All other trademarks are the sole property of their respective owners.

Part Number 4401661 Rev. C 06/2010

# Contents

	Preface xiii
	Safety information
	About the product
	Purpose of this guide
	Audience
	Assumptions
	How to use this guide
	How to obtain support xvii
Chapter 1	Instrument and Software Description
	System description
	Instrument description
	Instrument interior components
	Instrument parts and functions
	Theory of operation
	Preparing samples
	Preparing the instrument
	During a run
	Results
	Normalization
	Overview of the normalization feature7
	When to use the normalization feature 7
	Materials for routine operation
	External barcode scanner
	Uninterruptible Power Supply (UPS) 8
	Instrument reagents and consumables 9
	Anode buffer container (ABC)
	Cathode buffer container (CBC) 9
	Polymers
	Conditioning reagent
	Hi-Di <sup>™</sup> Formamide
	Capillary arrays
	Overview of the 3500 Series Data Collection Software
	About the software
	Parts of the software

	Use the software without an instrument	19
Chapter 2	Start the System	21
	Workflow	21
	Start the instrument	
	Start the computer	
	Log on to Windows	
	Launch the application	
	Log In	
	Check system status in the Dashboard	
	Dashboard, a quick glance	
	Check maintenance notifications	
	Check consumable status	
	Check buffer fill levels	31
	Replenish consumables	31
	Set preferences	
	Overview	32
	System preferences	32
	User preferences	33
	Set general preferences	33
	Set sequencing preferences	36
Chapter 3	Set Up and Run	41
	Workflow	
	Prepare the instrument	
	Create a plate	
	Assign plate contents	
	Name samples and assign sample types in the plate view	
	Assign assay, file name convention, and results group in the Plate View	
	Print the plate layout	
	Prepare and load sample plates	
	Prepare sample plates	
	Prepare the plate assembly	
	Load the plate in the instrument	
	Check instrument status	
	Link the plate	
	Quick Start a run	
	Load plates for run and create the injection list	
	Review and modify the injection list in Preview Run	
	Start the run	
	Monitor the run	61
	Check sequence or sample quality and specify re-injections	62

Specify re-injections       65         Review completed injections in Review Results       69         Start and stop a run       69         More features in Assign Plate Contents       70         Use the Plate View       70         Use the Table View       70         Use the Table View       71         Sort and customize tables       72         Add assays, file name conventions, and results groups to a plate       73         Create a plate for importing       73         Edit a plate       74         Import and export a plate       75         Specify the default plate type for the Open Plate dialog box       76         Save electronic version of reports       76         More features in Load Plate for Run       76         Link a plate from the recent plates or recent runs tab       76         More features in Monitor Run       77         Review traces       90         Review Sequencing Results       79         Review sequence quality       61         Review sequencing Results screen       80         Access the View Fagment/HID Analysis results       86         View, print, and save (.pdf) trace quality reports       85         Export sequencing results       87 <tr< th=""><th></th><th>Check sequence or sample quality</th><th> 62</th></tr<>		Check sequence or sample quality	62
Review completed injections in Review Results       69         Start and stop a run       69         More features in Assign Plate Contents       70         Use the Plate View       71         Sort and customize tables       72         Add assays, file name conventions, and results groups to a plate       73         Create a plate for importing       73         Edit a plate       74         Import and export a plate       75         Specify the default plate type for the Open Plate dialog box       76         Save electronic version of reports       76         More features in Load Plate for Run       76         More features in Monitor Run       77         Review Results       79         Review Sequencing Results       80         Access the View Sequencing Results screen       80         Review traces       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (pdf) trace quality reports       85         Review sample quality       89         Review mormalized data       90         Review in one screent Plate screen       86         Noter starters in Review Results       97         Review t			
Start and stop a run       69         More features in Assign Plate Contents       70         Use the Plate View       70         Use the Table View       70         Use the Table View       71         Sort and customize tables       72         Add assays, file name conventions, and results groups to a plate       73         Create a plate for importing       73         Edit a plate       74         Import and export a plate       75         Specify the default plate type for the Open Plate dialog box       76         Save electronic version of reports       76         More features in Load Plate for Run       76         Link a plate from the recent plates or recent runs tab       76         More features in Monitor Run       77         Review Results       79         Review Sequencing Results screen       80         Access the View Sequencing Results screen       80         Access the View Sequencing Results       87         Review sequence quality       81         Review sequence quality       84         Specify re-injections       85         View, print, and save (pdf) trace quality reports       86         Access the View Fragment/HID Results screen       88			
More features in Assign Plate Contents       70         Use the Plate View       70         Use the Table View       71         Sort and customize tables       72         Add assays, file name conventions, and results groups to a plate       73         Create a plate for importing       73         Edit a plate       74         Import and export a plate       75         Create a plate template       75         Specify the default plate type for the Open Plate dialog box       76         Save electronic version of reports       76         More features in Load Plate for Run       76         Link a plate from the recent plates or recent runs tab       76         More features in Monitor Run       77         Review the Instrument Run views       77         Review Sequencing Results       80         Access the View Sequencing Results screen       80         Review sequence quality       81         Review traces       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (pdf) trace quality reports       85         Review sample quality       89         Review sample quality       89         Revie			
Use the Plate View			
Use the Table View		-	
Sort and customize tables       72         Add assays, file name conventions, and results groups to a plate       73         Create a plate for importing       73         Edit a plate       74         Import and export a plate       75         Create a plate template       75         Specify the default plate type for the Open Plate dialog box       76         Save electronic version of reports       76         More features in Load Plate for Run       76         Link a plate from the recent plates or recent runs tab       76         More features in Load Plate for Run       77         Review the Instrument Run views       77         Review the Instrument Run views       77         Review Sequencing Results       80         Access the View Sequencing Results screen       80         Access the View Sequencing Results       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (.pdf) trace quality reports       85         Export sequencing results       87         Review sample quality       89         Review sing equality       89         Review fragment/HID Analysis results screen       86         Review sizing <td< td=""><td></td><td></td><td></td></td<>			
Add assays, file name conventions, and results groups to a plate       73         Create a plate for importing       73         Edit a plate       74         Import and export a plate       75         Create a plate template       75         Specify the default plate type for the Open Plate dialog box       76         Save electronic version of reports       76         More features in Load Plate for Run       76         Link a plate from the recent plates or recent runs tab       76         More features in Monitor Run       77         Review the Instrument Run views       77         Review sequencing Results       80         Access the View Sequencing Results screen       80         Review sequence quality       81         Review traces       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (pdf) trace quality reports       85         Export sequencing results       86         Access the View Fragment/HID Results screen       88         Review races       87         Export sequencing results       87         Review normalized data       90         Review protes       91         Review			
Create a plate for importing       73         Edit a plate       74         Import and export a plate       75         Create a plate template       75         Specify the default plate type for the Open Plate dialog box       76         Save electronic version of reports       76         More features in Load Plate for Run       76         Link a plate from the recent plates or recent runs tab       76         More features in Monitor Run       77         Review the Instrument Run views       77         Review the Instrument Run views       77         Review Sequencing Results       80         Access the View Sequencing Results screen       80         Review sequence quality       81         Review traces       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (.pdf) trace quality reports       85         Export sequencing results       87         Review sample quality       89         Review mormalized data       90         Review normalized data       90         Review sizing       94         Specify re-injections       95         View, print, and save (.pdf) sample quality reports <td></td> <td></td> <td></td>			
Edit a plate       74         Import and export a plate       75         Create a plate template       75         Specify the default plate type for the Open Plate dialog box       76         Save electronic version of reports       76         More features in Load Plate for Run       76         Link a plate from the recent plates or recent runs tab       76         More features in Monitor Run       77         Review the Instrument Run views       77         Chapter 4       Review Results       79         Review Sequencing Results       80         Access the View Sequencing Results screen       80         Access the View Sequencing Results screen       81         Review traces       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (,pdf) trace quality reports       85         Review sample quality       89         Review sample quality       89         Review sample quality       89         Review sample quality       89         Review protes       91         Review sing       91         Review sing       94         Specify re-injections       95 <t< td=""><td></td><td></td><td></td></t<>			
Import and export a plate       75         Create a plate template       75         Specify the default plate type for the Open Plate dialog box       76         Save electronic version of reports       76         More features in Load Plate for Run       76         Link a plate from the recent plates or recent runs tab       76         More features in Monitor Run       77         Review Results       77         Review Sequencing Results       80         Access the View Sequencing Results screen       80         Review traces       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (.pdf) trace quality reports       85         Export sequencing results       88         Access the View Fragment/HID Analysis results screen       88         Review sample quality       89         Review plots       91         Review plots       91         Review sizing       94         Specify re-injections       95         View, print, and save (.pdf) sample quality reports       85         Review sample quality       89         Review sample quality       89         Review sample quality       89<			
Create a plate template       75         Specify the default plate type for the Open Plate dialog box       76         Save electronic version of reports       76         More features in Load Plate for Run       76         Link a plate from the recent plates or recent runs tab       76         More features in Monitor Run       77         Review the Instrument Run views       77         Review Sequencing Results       80         Access the View Sequencing Results screen       80         Review sequence quality       81         Review traces       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (pdf) trace quality reports       85         Export sequencing results       88         Access the View Fragment/HID Results screen       88         Review sample quality       89         Review sizing       91         Review sizing       94         Specify re-injections       95         View, print, and save (pdf) sample quality reports       95         Review sample quality       89         Review sample quality       89         Review sample quality       89         Review plots <t< td=""><td></td><td>•</td><td></td></t<>		•	
Specify the default plate type for the Open Plate dialog box       76         Save electronic version of reports       76         More features in Load Plate for Run       76         Link a plate from the recent plates or recent runs tab       76         More features in Monitor Run       77         Review the Instrument Run views       77         Review Results       79         Review Sequencing Results       80         Access the View Sequencing Results screen       80         Review sequence quality       81         Review traces       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (pdf) trace quality reports       85         Export sequencing results       88         Access the View Fragment/HID Analysis results       88         Access the View Fragment/HID Results screen       88         Review sample quality       89         Review normalized data       90         Review sizing       94         Specify re-injections       95         View, print, and save (pdf) sample quality reports       95         Review sizing       94         Specify re-injections       95         View, prin			
Save electronic version of reports       76         More features in Load Plate for Run       76         Link a plate from the recent plates or recent runs tab       76         More features in Monitor Run       77         Review the Instrument Run views       77         Review Sequencing Results       80         Access the View Sequencing Results screen       80         Review sequence quality       81         Review traces       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (.pdf) trace quality reports       85         Export sequencing results       87         Review sample quality       88         Access the View Fragment/HID Analysis results       88         Review sizing       91         Review sizing       91         Review sizing       94         Specify re-injections       95         View, print, and save (.pdf) sample quality reports       95         Export secure of the sample quality       95         Review sizing       95         Review plots       95         Nore features in Review Results       97         Use Rename       97         S			
More features in Load Plate for Run       76         Link a plate from the recent plates or recent runs tab       76         More features in Monitor Run       77         Review the Instrument Run views       77         Chapter 4       Review Results       79         Review Sequencing Results       80         Access the View Sequencing Results screen       80         Review sequence quality       81         Review traces       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (.pdf) trace quality reports       85         Export sequencing results       88         Access the View Fragment/HID Analysis results screen       88         Review sample quality       89         Review sizing       91         Review sizing       94         Specify re-injections       95         View, print, and save (.pdf) sample quality reports       95         Export scip results       96         More features in Review Results       97         Use Rename       97         Sort       97			
Link a plate from the recent plates or recent runs tab       76         More features in Monitor Run       77         Review the Instrument Run views       77         Chapter 4       Review Results       79         Review Sequencing Results       80         Access the View Sequencing Results screen       80         Access the View Sequencing Results screen       80         Review sequence quality       81         Review races       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (.pdf) trace quality reports       85         Export sequencing results       87         Review Fragment/HID Analysis results       88         Access the View Fragment/HID Results screen       88         Review sample quality       89         Review wormalized data       90         Review sizing       94         Specify re-injections       95         View, print, and save (.pdf) sample quality reports       95         Keview pint, and save (.pdf) sample quality reports       95         Nore features in Review Results       97         Use Rename       97         Sort       97			
More features in Monitor Run       77         Review the Instrument Run views       77         Chapter 4       Review Results       79         Review Sequencing Results       80         Access the View Sequencing Results screen       80         Review sequence quality       81         Review traces       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (.pdf) trace quality reports       85         Export sequencing results       87         Review Fragment/HID Analysis results       88         Access the View Fragment/HID Results screen       88         Review somple quality       89         Review normalized data       90         Review sizing       94         Specify re-injections       95         View, print, and save (.pdf) sample quality reports       95         Nerview ports       95         More features in Review Results       97         Use Rename       97         Sort       97			
Review the Instrument Run views       77         Chapter 4       Review Results       79         Review Sequencing Results       80         Access the View Sequencing Results screen       80         Review sequence quality       81         Review traces       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (.pdf) trace quality reports       85         Export sequencing results       87         Review Fragment/HID Analysis results       88         Access the View Fragment/HID Results screen       88         Review sample quality       89         Review plots       91         Review sizing       94         Specify re-injections       95         View, print, and save (.pdf) sample quality reports       95         More features in Review Results       97         Use Rename       97         Sort       97			
Chapter 4       Review Results       79         Review Sequencing Results       80         Access the View Sequencing Results screen       80         Review sequence quality       81         Review traces       82         Understand Quality Values (QVs)       84         Specify re-injections       85         View, print, and save (.pdf) trace quality reports       85         Export sequencing results       87         Review Fragment/HID Analysis results       88         Access the View Fragment/HID Results screen       88         Review sample quality       89         Review plots       91         Review plots       91         Review sizing       94         Specify re-injections       95         View, print, and save (.pdf) sample quality reports       95         View, print, and save (.pdf) sample quality reports       95         View, print, and save (.pdf) sample quality reports       95         View, print, and save (.pdf) sample quality reports       95         Kiew print, and save (.pdf) sample quality reports       95         Kiew print, and save (.pdf) sample quality reports       95         Kiew print, and save (.pdf) sample quality reports       95         Kiew print, and			
Review Sequencing Results80Access the View Sequencing Results screen80Review sequence quality81Review traces82Understand Quality Values (QVs)84Specify re-injections85View, print, and save (.pdf) trace quality reports85Export sequencing results87Review Fragment/HID Analysis results88Access the View Fragment/HID Results screen88Review sample quality89Review plots91Review plots91Review sizing94Specify re-injections95View, print, and save (.pdf) sample quality reports95Export sizing results96More features in Review Results97Use Rename97Sort97			
Access the View Sequencing Results screen80Review sequence quality81Review traces82Understand Quality Values (QVs)84Specify re-injections85View, print, and save (.pdf) trace quality reports85Export sequencing results87Review Fragment/HID Analysis results88Access the View Fragment/HID Results screen88Review sample quality89Review plots91Review sizing94Specify re-injections95View, print, and save (.pdf) sample quality reports95Export sizing results96More features in Review Results97Use Rename97Sort97	Chapter 4	Review Results	79
Access the View Sequencing Results screen80Review sequence quality81Review traces82Understand Quality Values (QVs)84Specify re-injections85View, print, and save (.pdf) trace quality reports85Export sequencing results87Review Fragment/HID Analysis results88Access the View Fragment/HID Results screen88Review sample quality89Review plots91Review sizing94Specify re-injections95View, print, and save (.pdf) sample quality reports95Export sizing results97Use Rename97Sort97		Review Sequencing Results	80
Review traces82Understand Quality Values (QVs)84Specify re-injections85View, print, and save (.pdf) trace quality reports85Export sequencing results87Review Fragment/HID Analysis results88Access the View Fragment/HID Results screen88Review sample quality89Review normalized data90Review plots91Review sizing94Specify re-injections95View, print, and save (.pdf) sample quality reports95Export sizing results96More features in Review Results97Use Rename97Sort97			
Review traces82Understand Quality Values (QVs)84Specify re-injections85View, print, and save (.pdf) trace quality reports85Export sequencing results87Review Fragment/HID Analysis results88Access the View Fragment/HID Results screen88Review sample quality89Review normalized data90Review plots91Review sizing94Specify re-injections95View, print, and save (.pdf) sample quality reports95Export sizing results96More features in Review Results97Use Rename97Sort97			
Specify re-injections85View, print, and save (.pdf) trace quality reports85Export sequencing results87Review Fragment/HID Analysis results88Access the View Fragment/HID Results screen88Review sample quality89Review normalized data90Review plots91Review sizing94Specify re-injections95View, print, and save (.pdf) sample quality reports95Export sizing results96More features in Review Results97Use Rename97Sort97			
Specify re-injections85View, print, and save (.pdf) trace quality reports85Export sequencing results87Review Fragment/HID Analysis results88Access the View Fragment/HID Results screen88Review sample quality89Review normalized data90Review plots91Review sizing94Specify re-injections95View, print, and save (.pdf) sample quality reports95Export sizing results96More features in Review Results97Use Rename97Sort97		Understand Quality Values (QVs)	84
Export sequencing results87Review Fragment/HID Analysis results88Access the View Fragment/HID Results screen88Review sample quality89Review normalized data90Review plots91Review sizing94Specify re-injections95View, print, and save (.pdf) sample quality reports95Export sizing results96More features in Review Results97Use Rename97Sort97			
Export sequencing results87Review Fragment/HID Analysis results88Access the View Fragment/HID Results screen88Review sample quality89Review normalized data90Review plots91Review sizing94Specify re-injections95View, print, and save (.pdf) sample quality reports95Export sizing results96More features in Review Results97Use Rename97Sort97		View, print, and save (.pdf) trace quality reports	85
Access the View Fragment/HID Results screen88Review sample quality89Review normalized data90Review plots91Review sizing94Specify re-injections95View, print, and save (.pdf) sample quality reports95Export sizing results96More features in Review Results97Use Rename97Sort97			
Review sample quality89Review normalized data90Review plots91Review sizing94Specify re-injections95View, print, and save (.pdf) sample quality reports95Export sizing results96More features in Review Results97Use Rename97Sort97		Review Fragment/HID Analysis results	88
Review sample quality89Review normalized data90Review plots91Review sizing94Specify re-injections95View, print, and save (.pdf) sample quality reports95Export sizing results96More features in Review Results97Use Rename97Sort97		Access the View Fragment/HID Results screen	88
Review normalized data90Review plots91Review sizing94Specify re-injections95View, print, and save (.pdf) sample quality reports95Export sizing results96More features in Review Results97Use Rename97Sort97		Review sample quality	89
Review sizing94Specify re-injections95View, print, and save (.pdf) sample quality reports95Export sizing results96More features in Review Results97Use Rename97Sort97			
Specify re-injections       95         View, print, and save (.pdf) sample quality reports       95         Export sizing results       96         More features in Review Results       97         Use Rename       97         Sort       97		Review plots	91
Specify re-injections       95         View, print, and save (.pdf) sample quality reports       95         Export sizing results       96         More features in Review Results       97         Use Rename       97         Sort       97		Review sizing	94
Export sizing results		-	
Export sizing results		View, print, and save (.pdf) sample quality reports	95
More features in Review Results       97         Use Rename       97         Sort       97			
Use Rename			
Sort			
······································		Modify sequence, fragment analysis, or HID data	
		More features in Review Results	97

Chapter 5	Calibrate and Check Performance	. 99
	Section 1 Calibration	99
	Spatial calibration	99
	When to perform a spatial calibration	99
	Perform a spatial calibration	99
	Evaluate the spatial calibration profile	101
	Example spatial profiles	101
	Export spatial calibration results	101
	View and print a spatial calibration report	102
	Save historical calibration reports (.pdf) for record keeping	102
	Spectral calibration	103
	When to perform a spectral calibration	103
	Prepare for the spectral calibration	104
	Perform a spectral calibration	106
	Evaluate the spectral calibration data	110
	What you see during a spectral calibration	112
	Example spectral calibration data	114
	Export spectral calibration results	115
	View and print a spectral calibration report	116
	Save historical calibration reports (.pdf) for record keeping	116
	View the spectral calibration history	117
	Section 2 Performance check	. 119
	Run the sequencing install standard performance check	119
	Prepare for the sequencing install standard performance check	119
	Run the sequencing install standard performance check	122
	What you see during a run	124
	How the software determines passing and failing capillaries for the spectral calib 125	oration
	How the software determines passing and failing capillaries for the sequencing	
	performance check	
	Example sequencing install standard results	
	View previously run sequencing install standards	
	View and print a sequencing install standard report	
	Save historical performance check reports (.pdf) for record keeping	
	Run the fragment analysis or HID Install standard performance check	
	When to perform	
	Prepare for the fragment or HID install standard performance check	
	Run the fragment analysis or HID install standard performance check	
	What you see during a run	
	How the software determines passing and failing capillaries for the fragment/HI	
	performance check	

	Evaluate fragment/HID install standard data	135
	Example fragment install standard results	137
	Example HID install standard results	137
	View previously run install standards	137
	View and print a fragment or HID install standard detail report	137
	Save historical performance check reports (.pdf) for record keeping	138
Chapter 6	Manage Library Resources	. 139
	Overview of libraries	139
	General library procedures	140
	Access libraries	
	Create a new entry from a factory-provided, template, or locked entry	
	Delete a library entry	
	Edit a library entry	
	Import and export a library entry	
	View audit and e-signature histories for library entries	
	Sort, filter, and search library entries	
	Customize a library table	
	Plates library	
	Plate overview	
	Create a new plate	144
	Assays library	
	Assay overview	
	Create a new assay	
	File name conventions library	
	File name convention overview	
	Create a new file name convention	151
	Result group library	
	Results group overview	
	Create a new results group	
	Results group example 1: store files by plate name	
	Results group example 2: store one allelic ladder per run folder (8-capillary instruments)	
	Results group example 3: store re-injections in separate folders	162
	Instrument protocol library	
	Instrument protocol overview	165
	Create a new instrument protocol	
	Dye sets library	168
	Dye set overview	168
	Create a new dye set	
	Size standards library	
	Size standard overview	
	Normalization size standards provided	171

Create a new size standard	172
Modify a factory-provided normalization size standard	173
Basecalling protocols library (primary analysis – sequencing)	174
Basecalling protocol overview	174
Create a new basecalling protocol	174
Sizecalling protocols library (primary analysis – fragment)	179
Sizecalling protocol overview	179
Create a new sizecalling protocol	179
QC protocols library (primary analysis – HID)	184
QC protocol overview	184
Create a new QC protocol	184
Sequencing analysis protocols library (secondary analysis)	189
Sequencing analysis protocol overview	189
Create a new sequencing analysis protocol	189
MicroSeq <sup>®</sup> ID protocols library (secondary analysis)	
MicroSeq <sup>®</sup> ID analysis protocol overview	
Create a new MicroSeq <sup>®</sup> ID analysis protocol	
Fragment analysis protocols library (secondary analysis)	
Fragment analysis protocol overview	
Create a new fragment analysis protocol	
HID analysis protocols library (secondary analysis)	
HID analysis protocol overview	
Create a new HID analysis protocol	195

## Chapter 7 Use Security, Audit, and E-Sig Functions (SAE Module) ... 197

Section 1 Administrators 197
Administrators overview of system security, auditing, and electronic signature 197
Configure the security system 198
Access the Security screen and enable or disable security
Set account setup and security policies
Set up messaging notifications 200
Manage user accounts
Create or edit a user account
Determine the name of the logged-in user
Create or edit a user role
View and print a user report
Manage auditing
Access the Audit screen and enable or disable auditing
Select objects to audit
Create audit reason settings 208
Generate audit reports
Display audit histories

	Review the object audit history	210
	Review the system configuration history	211
	Review the action log	212
	View and print audit reports	213
	Archive, purge, and restore audit records	214
	Export audit records	215
	Manage electronic signature	216
	Access the E-Signature Settings screen and enable or disable e-sig	216
	Select the actions that allow signature	217
	Generate e-signature reports	221
	Display e-signature records	221
	View and print e-signature reports	222
	Export e-sig records	223
	Export and import user accounts, security, audit, and electronic signature settings	224
	Section 2 Users	. 225
	Users overview of System Security, Audit Trail and E-Signature	225
	Security	
	Audit	227
	Electronic signature	227
Ole and an O		000
Chapter 8	Maintain the Instrument	229
	Maintenance schedule	229
	Review maintenance notifications	229
	Daily instrument maintenance tasks	230
	Weekly instrument maintenance tasks	231
	Monthly instrument maintenance tasks	231
	Quarterly maintenance tasks	231
	Annual planned maintenance tasks	232
	As-needed instrument maintenance tasks	232
	Use the maintenance calendar	232
	View the calendar	232
	Default calendar entries	233
	Create calendar entries	233
	Review the Maintenance Notifications Log	234
	Instrument operational procedures	236
	Check consumables on the Dashboard	236
	Change the anode buffer container (ABC)	237
	Change the cathode buffer container (CBC)	238
	Check stored capillary arrays	240
		0/1
	Flush the water trap (pump trap)	241
	Flush the water trap (pump trap)          Routine instrument cleaning	

	Use the Maintenance Wizards to perform operations	. 244
	About Maintenance Wizards	. 244
	Replenish polymer	. 245
	Change polymer type	. 247
	Partially used polymer	. 249
	Wash the pump chamber and channels	. 249
	Use the conditioning reagent	. 250
	Fill capillary array with fresh polymer	. 251
	Remove bubbles from the polymer pump	. 251
	To change the capillary array	. 252
	To shutdown the instrument	. 253
	Computer maintenance	. 254
	Uninstall the software	. 254
	Archive, purge, and restore data	. 254
	Monitor disk space	. 256
	Review the Maintenance Notifications Log	. 257
	Service Log	. 258
Appendix A	Application Reagents and Run Modules	259
	Sequencing analysis reagents	. 259
	Fragment and HID analysis reagents	. 260
	Sequencing analysis dye sets for all applications	. 261
	Fragment analysis dye sets for all applications	. 261
	HID analysis dye sets	. 262
	Run modules	. 263
	Capillary array and polymer (sequencing analysis run modules)	. 263
	Capillary array and polymer (fragment and HID analysis run modules)	. 264
A		007
Appendix B	Secondary Analysis: Sequencing	267
	Perform secondary analysis on sequencing experiments	. 267
	Auto-analyze projects in the sequencing analysis software	. 267
	Set up an auto-analysis project in SeqScape®	. 268
	Set up a SeqScape plate in the 3500 Series Data Collection Software	. 270
	Auto-analysis with MicroSeq® ID	. 277
Appendix C	Secondary Analysis: Fragment	279
	Perform secondary analysis on fragment experiments	
	Auto-analyze projects in the fragment analysis software	
	Set up an auto-analysis project in GeneMapper®	
	Set up a GeneMapper plate in the 3500 Series Data Collection Software	
	Auto-Analysis with GeneMapper® ID-X	. 289

Appendix D	Remote Auto-Analysis Setup	291
	Remote auto-analysis configuration	291
	Remote auto-analysis installation	
	Create a shared folder	
	Set up the 3500 Series Data Collection Software v1.0	297
Appendix E	Troubleshoot	299
	Instrument troubleshooting	299
	Spatial calibration troubleshooting	
	Spectral calibration troubleshooting	
	Sequencing install standard troubleshooting	
	Fragment/HID install standard troubleshooting	
	Anode buffer container troubleshooting	303
	Cathode buffer container troubleshooting	304
	RFID troubleshooting	304
	Link a plate troubleshooting	304
	How to search and use the log files	305
	View instrument sensor details	305
	Data/electropherogram troubleshooting	306
	Dashboard troubleshooting	309
	Load plate troubleshooting	309
	Monitor run troubleshooting	311
	Review results troubleshooting	311
	Review error message details	312
	Audit troubleshooting	312
	Electronic signature troubleshooting	312
	Manual commands troubleshooting	313
	Miscellaneous	313
	Reset the instrument	314
		o / F
Appendix F	Safety	315
	Instrumentation safety	315
	Symbols on instruments	315
	Safety labels on instruments	317
	General instrument safety	317
	Physical hazard safety	318
	Electrical safety	319
	Laser safety	320
	Bar code scanner laser safety	321
	Workstation safety	321
	Safety and electromagnetic compatibility (EMC) standards	322

Chemical safety
General chemical safety 328
MSDSs
Chemical waste safety
Biological hazard safety
Safety alerts
Chemical alerts
Instrumentation alerts
Documentation
Related documentation
Obtaining information from the Help system
Send us your comments
Index

# Preface

## Safety information

**Note:** For general safety information, see this Preface and Appendix F, "Safety" on page 315. When a hazard symbol and hazard type appear by a chemical name or instrument hazard, see Appendix F, "Safety" on page 315 for the complete alert on the chemical or instrument.

#### Safety alert words

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

**IMPORTANT!** – Indicates information that is necessary for proper instrument operation, accurate chemistry kit use, or safe use of a chemical.

**CAUTION!** – Indicates a potentially hazardous situation that, if not avoided, may result in minor or moderate injury. It may also be 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 IMPORTANTs, each safety alert word in an Applied Biosystems 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 Applied Biosystems instruments* (see "Safety symbols" on page 315).

**MSDSs** The MSDSs (Material Safety Data Sheets) for any chemicals supplied by Applied Biosystems or Ambion are available to you free 24 hours a day. For instructions on obtaining MSDSs, see "MSDSs" on page 329.

**IMPORTANT!** For the MSDSs of chemicals not distributed by Applied Biosystems or Ambion contact the chemical manufacturer.

# Safety labels on<br/>instrumentsThe following CAUTION, WARNING, and DANGER statements may be displayed<br/>on Applied Biosystems instruments in combination with the safety symbols<br/>described in the preceding section.

Hazard symbol	English	Français
	<b>CAUTION!</b> Hazardous chemicals. Read the Material Safety Data Sheets (MSDSs) before handling.	<b>ATTENTION!</b> Produits chimiques dangereux. Lire les fiches techniques de sûreté de matériels avant toute manipulation de produits.
	<b>CAUTION!</b> Hazardous waste. Refer to MSDS(s) and local regulations for handling and disposal.	<b>ATTENTION!</b> Déchets dangereux. Lire les fiches techniques de sûreté de matériels et la régulation locale associées à la manipulation et l'élimination des déchets.
K	<b>CAUTION!</b> Potential slipping hazard.	<b>ATTENTION!</b> Risque potentiel d'avoir un sol glissant.
	CAUTION! Hot surface.	ATTENTION! Surface brûlante.
$\wedge$	DANGER! High voltage.	DANGER! Haute tension.
<u> </u>	<b>WARNING!</b> To reduce the chance of electrical shock, do not remove covers that require tool access. No user-serviceable parts are inside. Refer servicing to Applied Biosystems qualified service personnel.	<b>AVERTISSEMENT!</b> Pour éviter les risques d'électrocution, ne pas retirer les capots dont l'ouverture nécessite l'utilisation d'outils. L'instrument ne contient aucune pièce réparable par l'utilisateur. Toute intervention doit être effectuée par le personnel de service qualifié venant de chez Applied Biosystems.
*	<b>CAUTION!</b> Class 2(II) visible and/or invisible radiation present. Do not stare directly into the beam or view directly with optical instruments.	<b>ATTENTION!</b> Rayonnement visible ou invisible d'un faisceau. Ne pas regarder le faisceau directement ou au travers d'un instrument optique.
	<b>DANGER!</b> Class 3B (III) visible and/or invisible radiation present. Avoid exposure to beam.	<b>DANGER!</b> Rayonnement visible ou invisible d'un faisceau de Classe 3B (III) en cas d'ouverture. Evitez toute exposition au faisceau.
	CAUTION! Sharp object.	AVERTISSEMENT! Objet Sharp.

## About the product

The Applied Biosystems 3500/3500xL Genetic Analyzers is an automated 8 and/or 24 capillary instrument designed for a wide range of sequencing and fragment analysis applications.

IMPORTANT! For Research Use Only. Not for use in diagnostic procedures.

## Purpose of this guide

The *Applied Biosystems 3500/3500xL Genetic Analyzer User Guide* provides stepby-step instructions for preparing and analyzing a sample. It is designed to help you learn how to use the instrument.

**CAUTION!** The protection provided by the equipment may be impaired if the instrument is operated outside the environment and use specifications, the user provides inadequate maintenance, or the equipment is used in a manner not specified by the manufacturer (Applied Biosystems).

## Audience

This user guide is written for principle investigators and laboratory staff who are planning to operate and maintain the Applied Biosystems 3500/3500xL Genetic Analyzers.

## Assumptions

The *Applied Biosystems 3500/3500xL Genetic Analyzer User Guide* assumes that your 3500 or 3500xL analyzer has been installed by an Applied Biosystems service representative.

This guide also assumes you have the following background:

- Familiarity with Microsoft<sup>®</sup> Windows Vista<sup>®</sup> operating system.
- Knowledge of general techniques for handling DNA samples and preparing them for electrophoresis.
- A general understanding of hard drives, data storage, file transfers, and copying and pasting.

## How to use this guide

EPT

FR

NIC

NT

Pe

QV

MicroSeq<sup>®</sup> kit

(or other product name)

Text conventions	This guide uses the following conve	entions:					
	• <b>Bold</b> text indicates user action. For example:						
	Type <b>0</b> , then press <b>Enter</b> for each of the remaining fields.						
	<ul> <li><i>Italic</i> text indicates new or important words and is also used for emphasis.</li> <li>For example:</li> </ul>						
	Before changing reagents, <i>always</i> determine what chemicals have been used in the instrument.						
	<ul> <li>A right arrow symbol ( ) separates successive commands you select from a drop-down or shortcut menu. For example:</li> </ul>						
	Select File > Open > Spot Set.						
	Right-click the sample row, then select <b>View Filter &gt; View All Runs</b> .						
User attention words							
	<b>Note:</b> Provides information that may be of interest or help but is not critical to the use of the product.						
	<b>IMPORTANT!</b> Provides information that is necessary for proper instrument operation, accurate chemistry kit use, or safe use of a chemical.						
Table of Acronyms							
	Acronym	Definition					
	ABC	Anode Buffer Container					
	BDT	BigDye <sup>®</sup> Terminator Kit					
	BDX	BigDye <sup>®</sup> Xterminator <sup>™</sup>					
	Сар	Capillaries					
	CBC	Cathode Buffer Container					
	CV/Fitting Check Valve pouch attachment fitting						
	CV/Fitting Check Valve pouch attachment fitting						

Applied Biosystems 3500/3500xL Genetic Analyzer User Guide

Nucleotide Base Color (A, G, C, T)

**ElectroPhoresis Telemetry** 

**Factory Repeating** 

Nucleoid Type

**Quality Value** 

Probability of error

**Microbial Sequencing** 

Network Interface Card

Acronym	Definition			
GM	GeneMapper <sup>®</sup> Software			
GMIDx	GeneMapper <sup>®</sup> IDx			
POP™	Polymer (Brand name of the polymer)			
PPS	Power Protection System			
SAE	Security, Administration, Electronic signature			

## How to obtain support

For the latest services and support information for all locations, go to:

#### www.appliedbiosystems.com

At the Applied Biosystems web site, you can:

- Access worldwide telephone and fax numbers to contact Applied Biosystems Technical Support and Sales facilities.
- Search through frequently asked questions (FAQs).
- Submit a question directly to Technical Support.
- Order Applied Biosystems user documents, MSDSs, certificates of analysis, and other related documents.
- Download PDF documents.
- Obtain information about customer training.
- Download software updates and patches.

# How to obtain For detail more information 3500 Seri

For detailed information on preparing for installation, refer to the *Applied Biosystems* 3500 Series Genetic Analyzer Site Preparation Guide (4401689).

**Note:** The purpose of the Site Prep Guide is to help you prepare your site for installation of the 3500 or 3500xL analyzer. For specific details about your system, please refer to this user guide.

# Instrument and Software Description

## System description

The 3500 or 3500xL analyzer is shipped with the following system components:

- Capillary Electrophoresis instrument.
- 3500 (8-capillary) or 3500xL (24-capillary) array and POP<sup>™</sup> polymer
- DNA sequencing, or fragment analysis, reagents, and other consumables for system qualification.
- Dell<sup>®</sup> computer workstation with flat-screen monitor.
- Integrated software for instrument control, data collection, quality control, and basecalling or sizing of samples.

#### Instrument description

The Applied Biosystems 3500/3500xL Genetic Analyzers are fluorescence based DNA analysis instrument using capillary electrophoresis technology with 8- or 24-capillaries.

For detailed dimensions of the instrument, refer to the *Applied Biosystems 3500* Series Genetic Analyzer Site Preparation Guide (4401689).

**Note:** The purpose of the Site Prep Guide is to help you prepare your site for installation of the 3500 or 3500xL analyzer. For specific details about your system, please refer to this user guide.



### Instrument interior components

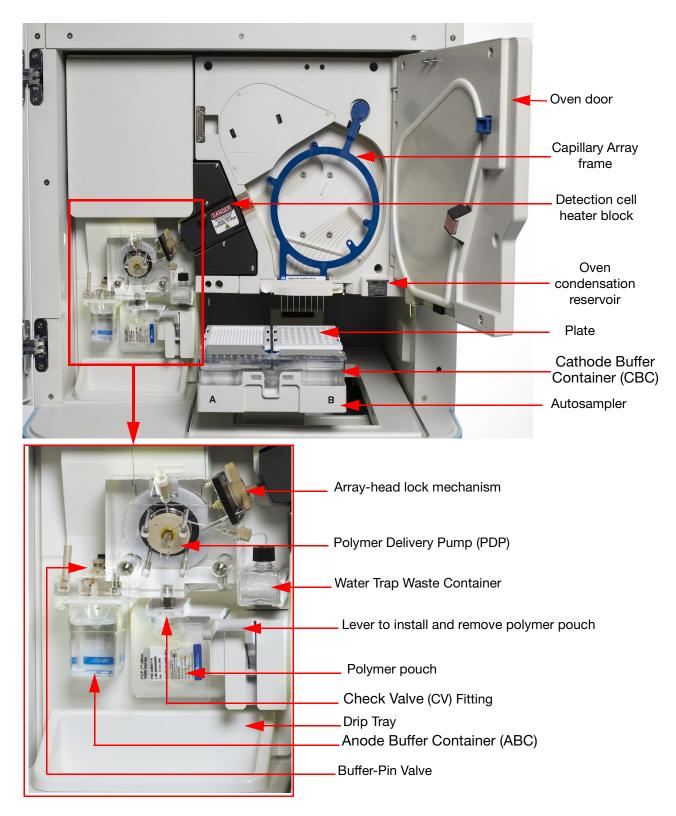


Figure 1 Instrument interior components

## Instrument parts and functions

#### Table 1 Instrument parts and functions

Part	Function
Autosampler	Holds the sample plates and Cathode Buffer Container (CBC) and moves to align the plates and CBC with the capillaries.
Oven	Maintains uniform capillary array temperature.
Oven condensation reservoir	Collects condensation from the oven.
Pump block	Includes the displacement pump chamber, polymer chambers, piston water seal, syringe fitting array attachment point (array port), the lower polymer block, and the CV/Fitting (Check Valve pouch attachment fitting).
Detection cell heater block	Holds the detection cell in place for laser detection and maintains the detection cell temperature of 50 °C.
Polymer Delivery Pump (PDP)	Pumps polymer into the array and allows for automated maintenance procedures.
Lower polymer block	Contains the buffer valve, anode electrode, buffer gasket, and holds the anode buffer container.
Radio Frequency Identification (RFID)	<ul> <li>RFID tags to read the following information for primary instrument consumables:</li> <li>Lot numbers</li> <li>Serial numbers</li> <li>Dates (expiration)</li> <li>Capacity (usage)</li> <li>The primary consumables are:</li> <li>Capillary Array</li> <li>Cathode Buffer Container (CBC)</li> <li>POP<sup>™</sup> Polymer</li> <li>Anode Buffer Container (ABC)</li> </ul>
Capillary Array	Enables the separation of the fluorescent-labeled DNA fragments by electrophoresis. It is a replaceable unit composed of 8 or 24 capillaries (50 cm and 36 cm length). Note: The 36 cm capillary is for HID applications, only.
Anode Buffer Container (ABC)	The Anode Buffer Container (ABC) contains 1× running buffer to support all electrophoresis applications on the instrument. It has a built-in overflow chamber to maintain constant fluid height.
Cathode Buffer Container (CBC)	The Cathode Buffer Container (CBC) contains 1× running buffer to support all electrophoresis applications on the instrument.
Polymer pouch	Supplies polymer to the Polymer Delivery Pump.
Conditioning reagent	The pouch is used for priming the polymer pump, washing the polymer pump between polymer type changes, and during instrument shut down. It has adequate volume for a one-time use.

## Theory of operation

The 3500 or 3500xL analyzer is a fluorescence-based DNA analysis system that uses proven capillary electrophoresis technology with 8- or 24-capillaries.

The 3500 or 3500xL analyzer is fully automated, from sample loading to primary data analysis, for sequencing, fragment analysis, and HID (human identification) analysis.

**Note:** In this document, primary analysis for sequencing is referred to as basecalling. Primary analysis for fragment or HID procedures is referred to as sizing.

#### **Preparing samples**

When DNA samples are prepared for sequencing, fragment analysis, or HID analysis on the 3500 or 3500xL analyzer, fluorescent dyes are attached to the DNA. For most applications, the sample is denatured so that only single-strand DNA is present.

#### Preparing the instrument

Two calibrations are required to prepare the instrument for sample runs:

- **Spatial calibration** Determines the position of the image from each capillary on the CCD array. For more information, refer to "Spatial calibration" on page 99.
- **Spectral calibration** Generates a matrix for each capillary that compensates for dye overlap and is used to convert the 20-color data into 4-, 5-, or 6-dye data. For more information, refer to "Spectral calibration" on page 103.

#### During a run

During a run, the system:

- Prepares the capillary by pumping fresh polymer solution under high pressure from the polymer delivery pump to the waste position in the Cathode Buffer Container (CBC).
- Electrokinetically injects the sample into the capillary using a low-voltage for a few seconds.
- Washes the capillary tips in the rinse position of the CBC, then returns the capillary to the buffer position of the CBC.

• Ramps the voltage up to a constant voltage.

A high electric field is created between the ground end of the Anode Buffer Container (ABC) and the negative voltage applied to the load header of the capillary array. This field pulls the negatively charged DNA through the separation polymer. The smaller fragments migrate faster than the larger fragments and reach the detector first.

To insure optimal separation and maintain denaturation of the DNA, the capillaries are thermally controlled in the oven and in the detection cell. The oven has a Peltier heat unit and fan-circulated air. The Peltier can heat and cool the oven to maintain sub-ambient temperatures, which are useful for non-denaturing applications such as SSCP (Single-strand conformation polymorphism).

- In the detection cell, the dyes attached to DNA are excited by a narrow beam of laser light. The laser light is directed into the plane of the capillaries from both the bottom and top. A small amount of laser light is absorbed by the dyes and emitted as longer wavelength light in all directions.
- Captures the fluorescent light on the instrument optics while blocking the laser light. The light passes through a transmission grating, which spreads the light out. The light is imaged onto a cooled, scientific-grade CCD array. For each capillary, 20 zones on the CCD are collected to provide20-color data for each capillary.
- Converts the 20-color data into multi-dye data for the entire run. For sequencing applications, 4 different dyes are used to determine the 4 bases A, G, C and T. For fragment and HID analysis applications, up to 6 dyes can be used in a single run for higher throughput.

#### Results

The software generates an electropherogram (intensity plot) for each dye based on the migration of DNA fragments over the run and generates primary analysis results:

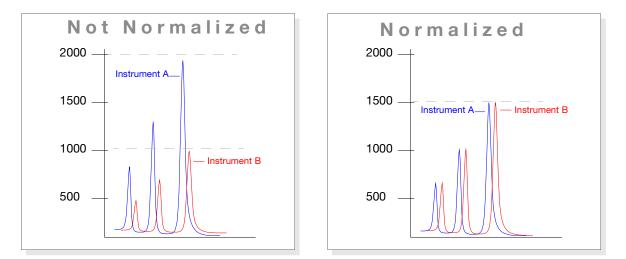
- For sequencing applications, the electropherogram is adjusted to compensate for slight mobility differences due to the dyes, then basecalling is performed and quality values are assigned.
- For fragment and HID analysis, the software uses the internal size standard to assign a fragment size and a sizing quality value to each peak.

If the autoanalysis functionality has been set up, the system transfers the sample data to a secondary analysis software application for further processing. Alternatively, you can manually transfer the sample data to a secondary analysis software application for further processing.

## Normalization

### Overview of the normalization feature

For fragment analysis and HID applications, the 3500 Series Data Collection Software includes a normalization feature for use with the GeneScan<sup>TM</sup> 600 LIZ<sup>®</sup> Size Standard v2.0 (GS600 LIZ v2). This feature attenuates signal variations associated with instrument, capillary array, sample salt load, and injection variability between capillaries and instruments. Normalization can be applied during primary analysis of the data.



# Figure 2 Comparison of fragment analysis results with and without the normalization feature

To use the normalization feature, prepare each sample with the GS600 LIZ v2 size standard, then specify the appropriate normalization size standard for file primary analysis. The GS600 LIZ v2 reagent can function as an internal standard for signal-height normalization as well as a size standard for peak sizing.

### When to use the normalization feature

The 3500 Series Data Collection Software provides three normalization size-standard definition files that you can specify for primary analysis of samples prepared with the GS600 LIZ v2 size standard:

- Fragment:
  - GS600LIZ+Normalization
  - GS600(60-600)LIZ+Normalization For applications that have primer peaks that obscure the 20 and 40-mer peaks of the GS600 LIZ v2 size standard.
- HID:
  - GS600(80-400)LIZ+Normalization

## Materials for routine operation

Contact your local Applied Biosystems service representative (or go to **www.appliedbiosystems.com**, then click **Products**) to order the materials for 3500 or 3500xL analyzer.

## External barcode scanner

An external barcode scanner can be used with the 3500 or 3500xL analyzer to scan the plate template.

Applied Biosystems recommends the Symbol LS 1203 handheld barcode scanner (shown), which connects to the instrument computer.

The scanner allows you to scan barcodes into any text box in the 3500 Series Data Collection Software.

For details on installation, use, decoding capabilities, and product specifications see the product documentation.

## Uninterruptible Power Supply (UPS)

Loss of power during a run can result in lost data. To address concerns with loss of power in the laboratories, Applied Biosystems recommends the use of an Instrumentation Power Protection System (IPPS) with the 3500 or 3500xL analyzer.

If your laboratory has a backup-power generator, a battery-powered backup Power Protection System (PPS) is required to provide power for the amount of time that it takes for your laboratory's backup power to begin generating power and stabilize.

Note: The instrument, computer, and monitor must all be connected to the PPS.

If your laboratory does not have a backup-power generator, a PPS can provide protection from power disruptions of a limited duration. For longer durations, optional battery cabinets can be added to the base PPS unit. A base unit PPS rated for 800W can provide over 20 minutes of backup protection, and over 2 hours when a single battery cabinet was added.

**Note:** Battery output can be affected by temperature and the age of battery so these backup times are not guaranteed.

## Instrument reagents and consumables

For application-specific reagents, consumables, and run modules, see Appendix A, Application Reagents and Run Modules.

#### Anode buffer container (ABC)

The ABC (PN 4393927) contains  $1 \times$  running buffer to support all electrophoresis applications on the 3500 or 3500xL analyzer.

The ABC is made in a ready to use, disposable, container with a radio frequency identification (RFID) tag incorporated into the label. It has a built-in overflow chamber to maintain constant fluid height.

For the following hazard(s), see the complete safety alert descriptions in "Specific chemical alerts" on page 333.

## WARNING! CHEMICAL HAZARD. Anode Buffer Container (ABC).

Store the ABC at 2 °C to 8 °C until ready to use. The sealed ABC is stable at this temperature until the expiration date shown on the label. Once the seal is peeled, the running buffer is stable at ambient temperature for up to 7 days. Ensure that the seal remains in place until just prior to use on the instrument.

To ensure optimal performance, the use of the ABC is limited to either 7 days after the first installation or 120 injections on a 3500 (8-capillary)/50 injections on a 3500xL (24-capillary), whichever comes first. When notified of the limit by the instrument software, you have to replace the ABC with a new one before you can proceed further.

For more details see the product insert included in the product package.

See "Change the anode buffer container (ABC)" on page 237 for instructions on how to change the ABC.

### Cathode buffer container (CBC)

The CBC (PN 4408256) contains  $1 \times$  running buffer to support all electrophoresis applications on the 3500 or 3500xL analyzer.

The CBC is made in a ready to use, disposable, container with a radio frequency identification (RFID) tag incorporated into the label. It has two separate sides:

- The side containing 24 holes provides the cathode buffer for electrophoresis.
- The side that contains 48 smaller holes provides the liquid for wash and waste functionality for rinsing the capillary tips and collecting wash waste between injections.

For the following hazard(s), see the complete safety alert descriptions in "Specific chemical alerts" on page 333.

#### WARNING! CHEMICAL HAZARD. Cathode Buffer Container (CBC).

Store the CBC at 2 °C to 8 °C until ready to use. The sealed CBC is stable at this temperature until the expiration date shown on the label. Once the seal is peeled, the running buffer is stable at ambient temperature for up to 7 days. Ensure that the seal remains in place until just prior to use on the instrument.

To ensure optimal performance, the use of the CBC is limited to either 7 days after the first installation or 120 injections on a 3500 (8-capillary)/50 injections on a 3500xL (24-capillary), whichever comes first. When notified of the limit by the instrument software, you have to replace the ABC with a new one before you can proceed further.

For more details see the product insert included in the product package.

See "Change the cathode buffer container (CBC)" on page 238 for instructions on how to change the CBC.

#### Polymers

The polymer for 3500 or 3500xL analyzer is available as a ready to use pouch with either POP-4<sup>TM</sup>, POP-6<sup>TM</sup> or POP-7<sup>TM</sup> polymer as the separation matrix.

**IMPORTANT!** Do not use a polymer pouch that has been installed on one type of instrument on another type of instrument. For example, if you install a new polymer pouch on a 3500 (8-capillary) instrument, do not use that polymer on a 3500xL (24-capillary) instrument.

The pouch has adequate polymer to support the stated number of samples (384 or 960) or injections and additional volume to handle a limited number of installations and setup related wizard operations. Incorporated into the label is a radio frequency identification (RFID) tag.

**Note:** The top part of the pouch fitment is sealed with a plastic film, which should be removed prior to direct installation on to the instrument.

For the following hazard(s), see the complete safety alert descriptions in "Specific chemical alerts" on page 333.

# **WARNING!** CHEMICAL HAZARD. POP- $4^{\text{TM}}$ , POP- $6^{\text{TM}}$ , and POP- $7^{\text{TM}}$ polymers.

Store the polymer at 2 °C to 8 °C until ready to use. The sealed polymer is stable at this temperature until the expiration date shown on the label.

For more details see the product insert included in the product package.

See "Change polymer type" on page 247 for instructions on how to change polymers.

**IMPORTANT!** If you remove a polymer pouch for storage, place a Pouch Cap (PN 4412619) onto the pouch, then place an empty pouch (or conditioning reagent) on the connector to prevent desiccation of any residual polymer on the connector. Follow the instructions in the wizard to ensure the proper operation of the pouch and the instrument.

#### Applications

- POP-6<sup>™</sup> and POP-7<sup>™</sup> polymers are recommended for sequencing and fragment analysis applications.
- POP-4<sup>TM</sup> polymer is recommended for HID/Forensic applications.

#### Table 2 Polymers used for all applications

				Pouch limits		
Polymer type	Part number	Instrument used	On-instrument life or whichever comes first <sup>‡</sup>	Cannot exceed <sup>§</sup>	User option to continue <sup>#</sup>	
POP-4 <sup>™</sup> (960) POP-6 <sup>™‡‡</sup> (960)	4393710 4393712	3500 (8-capillary)	Lower of 7 days or 960 samples or 120 Injections	Expiry date, Sample limit	7-day limit	
POP-7 <sup>™</sup> (960)	4393714	3500xL (24-capillary)	Lower of 7 days or 960 samples or 50 Injections	and/or Injections limit		
POP-4 <sup>™</sup> (384) POP-6 <sup>™ ‡‡</sup> (384)	4393715 4393717	3500 (8-capillary)	Lower of 7 days or 384 samples or 60 Injections	Expiry date, Sample limit	7-day limit	
POP-7 <sup>™</sup> (384)	4393708	3500xL (24-capillary)	Lower of 7 days or 384 samples or 20 Injections	and/or Injections limit		

The polymer pouch includes additional volume to accommodate a limited number of installation and wizard operations. However, if the number of wizard operations exceeds a certain limit, the number of remaining samples or injections will be reduced. For example, if you run the total bubble remove option in the bubble remove wizard more than four times, or run other wizards operations excessively, the number of remaining samples or injections will be reduced. Refer to the polymer gage on the dashboard for the up-to-date number of remaining samples or injections at any given point.

§ Replace the pouch before proceeding further.

# Applied Biosystems has verified the polymer for a maximum of 7 days on the instrument.

<sup>‡‡</sup> Ambient temperature must be in the range of 15 °C to 25 °C. Sustained use at higher temperatures may result in shorter read lengths than specified.

### Conditioning reagent

The conditioning reagent (PN 4393718) for 3500 or 3500xL analyzer is available as a ready to use pouch. It is used for priming the polymer pump, washing the polymer pump between polymer type changes, and during instrument shut down. It has adequate volume for a one-time use.

**Note:** Use of the conditioning reagent is dictated by the instrument wizards. Install the pouch when requested to do so by the wizard.

 $\Delta$  **CAUTION!** Expired pouches cannot be used on the instrument. Once installed on the instrument, the pouch is good for a one-time use, only.

For more details see the product insert included in the product package.

See "Use the conditioning reagent" on page 250 for instructions on how to use the conditioning reagent.

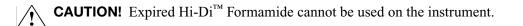
## Hi-Di<sup>™</sup> Formamide

Hi-Di<sup>™</sup> Formamide (pack of four) 5-ml tube (PN 4440753) is a highly deionized formamide, formulated with a stabilizer, ready for use as an injection solvent for all applications on the 3500 or 3500xL analyzer.

For the following hazard(s), see the complete safety alert descriptions in "Specific chemical alerts" on page 333.

#### WARNING! CHEMICAL HAZARD. Hi-Di<sup>™</sup> Formamide.

For more details see the product insert included in the product package.



#### Applications

The Hi-Di<sup>™</sup> Formamide is used for sequencing analysis, fragment analysis, and HID/Forensic applications. To determine the exact, and necessary, volume of formamide for each specific application, follow the provided protocols and product inserts.

Table 3 Hi-Di<sup>™</sup> Formamide used for all applications

Hi-Di <sup>™</sup> Formamide name	Instrument	Part number	On-instrument life and usage
Hi-Di <sup>™</sup> Formamide - 5-ml bottle (pack of four)	3500 (8-capillary) 3500xL (24-capillary)	4440753	24 hours

#### Capillary arrays

The capillary array for 3500 or 3500xL analyzer is installed on the instrument and ready to use.



**CAUTION!** SHARP The load-end of the capillary array has small but blunt ends and it could lead to piercing injury.

See "To change the capillary array" on page 252 for instructions on how to change the capillary array.

#### Applications

- The 36 cm capillary array is used for HID/Forensic applications.
- The 50 cm capillary array is used for sequencing and fragment analysis applications.

Table 4 Capillary arrays used for all applications
--

Capillary array name	Part number	Instrument used	On- instrument life	RFID-controlled limits	
			me	User option to continue <sup>‡</sup>	
8-Capillary, 36 cm	4404683	3500	160	Under user option to continue (160 injections	
8-Capillary, 50 cm	4404685	-	injections	and expiry date)	
24-Capillary, 36 cm	4404687	3500xL		Under user option to continue (160 injections	
24-Capillary, 50 cm	4404689	-	and expiry date)	and expiry date)	

‡ Applied Biosystems has verified the array for 160 injections.

# Overview of the 3500 Series Data Collection Software

## About the software

Application	Supports
Sequencing	<ul><li>Direct sequencing for mutation detection</li><li>Comparative sequencing with and without references</li><li>Microbial sequence identification</li></ul>
Fragment analysis	<ul> <li>Microsatellite</li> <li>AFLP<sup>®</sup> (amplified fragment length polymorphism)</li> <li>SNaPshot<sup>®</sup> kit</li> <li>LOH (loss of heterozygosity)</li> <li>MLPA<sup>®</sup> (Multiplex ligation-dependent probe amplification)</li> </ul>
HID	<ul><li>Forensic DNA casework</li><li>Databasing</li><li>Paternity testing</li></ul>

#### Table 5 3500 Series Data Collection Software applications supported

During a run, the software:

- Controls the instrument and generates sample data files:
  - Sequencing (.ab1)
  - Fragment analysis (.fsa)
  - HID (.hid)
- Performs primary analysis and reporting that evaluate the quality of the data:
  - Sequencing Basecalling and trimming
  - Fragment analysis and HID Peak detection and sizing
- (Optional) Performs secondary analysis (auto-analysis) with the following Applied Biosystems software applications:
  - Sequencing SeqScape<sup>®</sup> Software v2.7 (or later) or MicroSeq<sup>®</sup> ID Analysis Software v2.2 (or later)
  - Fragment analysis GeneMapper<sup>®</sup> Software v4.1 (or later)
  - HID GeneMapper<sup>®</sup> *ID-X* Software v1.2 (or later)

Note: You can also manually import sample data files in to the secondary analysis software applications above. Sample data files generated by the 3500 Series Data Collection Software are also compatible with Applied Biosystems Variant Reporter<sup>™</sup> Software (v1.1 or later) and Sequence Analysis (SeqA) Software (v5.4 or later).

### Parts of the software

**Dashboard** The first screen that is displayed when you start the 3500 Series Data Collection Software is the Dashboard (Figure 3).

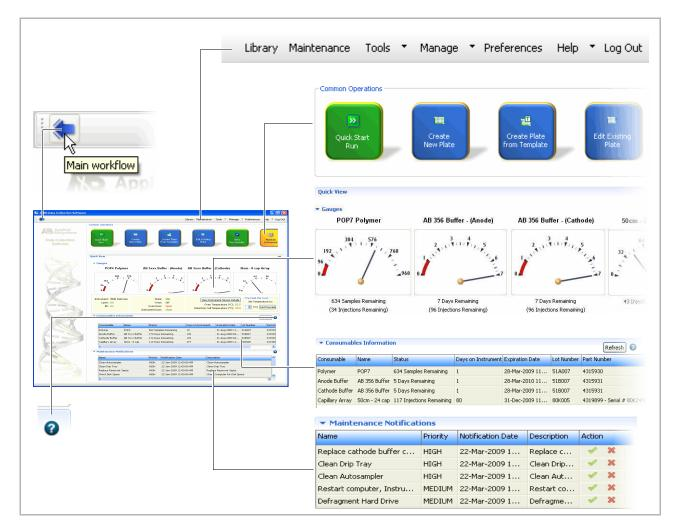


Figure 3 3500 Series Data Collection Software Dashboard

The Dashboard gives you quick access to the information and tasks you need to set up and run:

- Main workflow arrow Advances to the screens where you set up, load, and run plates, and view results.
- Menu bar Accesses all other parts of the software. The menu bar is displayed on all screens.
- **Common operations** Allows you to quick-start (load a plate that is set up), create or edit plates, view results, and access the Maintenance workflow.

- Quick view Displays gauges that show the remaining usage of consumables and gives the status of instrument conditions. Consumable usage is automatically tracked by the instrument by radio-frequencing identification (RFID) tags.
- **Consumables information** Gives details for the installed consumables and indicates in red if any consumable is about to expire based on RFID tags.
- Maintenance notifications Lists the scheduled maintenance tasks.
- Help icon ② Displays a help topic specific to a screen or an area of the screen.

For more information, see "Check system status in the Dashboard" on page 26.

**Main workflow** Click the main workflow arrow at the top left of the Dashboard to access the Main workflow.

The Main workflow contains the screens where you set up, load, and run plates, and view results.

The Main workflow navigation pane is designed as a task workflow. Each screen contains a button that you can click to advance to the next screen in the workflow.

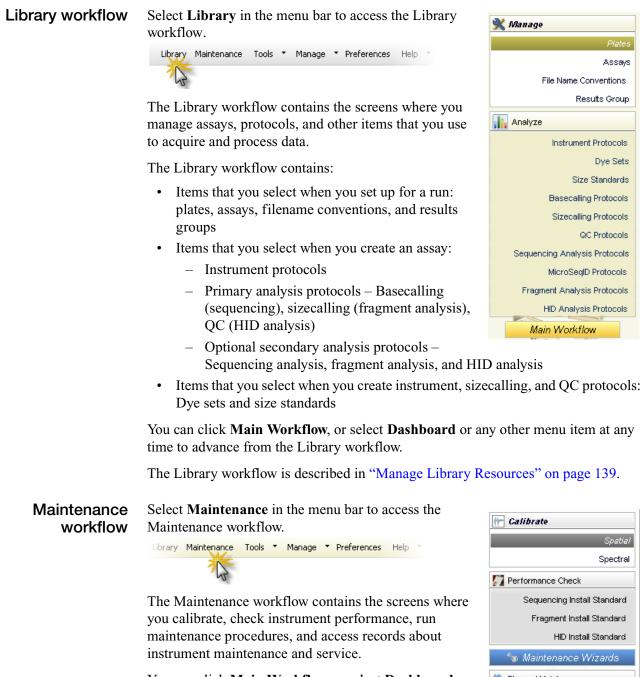
Select a task in the navigation pane to access each screen.

You can select **Dashboard** or any other menu item at any time to advance from the Main workflow.





The Main workflow is described in Chapter 3, "Set Up and Run" on page 41, and "Review Results" on page 79.



You can click Main Workflow or select Dashboard or any other menu item at any time to advance from the Maintenance workflow.

The Maintenance workflow is described in Chapter 8, "Maintain the Instrument" on page 229.

Sequencing Install Standard Fragment Install Standard HID Install Standard 物 Maintenance Wizards -Planned Maintenance Notifications Log Service Log Schedule Main Workflow

Tools menu	<ul> <li>Select Tools in the menu bar to access 3500 Series Data Collection Software tools.</li> <li>Tools provided are: <ul> <li>Security, Audit, and E-signature (if your system includes the SAE module)</li> <li>Change Password that allows you</li> <li>View Logs that provides reports of Manual Commands that you can be the SAE module is described in Chapp Functions (SAE Module)" on page 19</li> </ul> </li> </ul>	to cha of instr use to oter 7,	rument ru troublesh	ans. noot instrument	performance.
Manage menu	Select <b>Manage</b> in the menu bar to access archive, restore, and purge functions Archive, restore, and purge are described in Chapter 8, Maintain the Instrument.	Library M	laintenance	Tools • Manage Archive Restore Purge	Preferences Help
Preferences menu	<ul> <li>Select Preferences in the menu bar to access the parameters for which you can set defaults.</li> <li>Preferences allow you to set system and user defaults for settings st as the date format, sample data file storage location, export file formats for sequencing data, and a variety of sequencing-specific settings.</li> <li>System defaults apply to all users.</li> <li>User defaults apply to: <ul> <li>All users – If your system does n include the SAE module.</li> <li>Each logged-in user – If your system includes the SAE module.</li> </ul> </li> <li>Preferences are described in Chapter 2 "Start the System" on page 21.</li> </ul>	uch or ot	Le Prefer ype filter - Syster - Da - In: - Sc - Se - Se - Se - Se - Re - Ru	text not used	Preferences Help *  System Select one of the s system preferences enable you to do se
Help menu	Select <b>Help</b> in the menu bar to access 3500 Series Data Collection Software Help.	3	Preferences	Help Contents About 3500 Data C	ollection Software

The Help provides quick access to brief information about how to perform tasks on a screen. For details about tasks and other information, refer to the chapters in this user guide.

Navigate the Software

### From the Dashboard

To advance from the Dashboard to:

- Main workflow Click ().
- Other screens in the software Select items from the menu bar.

### From the Main workflow

To advance from the Main workflow to:

- Dashboard Click Dashboard.
- Other screens in the Main workflow Select items in the navigation pane.
- Other screens in the software Select items from the menu bar.

### From the Library or Maintenance workflows

To advance from the Maintenance or Library workflow to:

- Dashboard Click Dashboard.
- Other screens in the workflow Select items in the navigation pane.
- Main workflow Click Main Workflow in the navigation pane.
- Other screens in the software Select items from the menu bar.







### Use the software without an instrument

You can install the 3500 Series Data Collection Software on a computer that is not connected to an instrument. You can use this stand-alone version of the software to create plates, protocols, and other library items, and to review completed results.

**IMPORTANT!** Do not select instrument-related functions in the stand-alone version of the software.

# Start the System

## Workflow

## Start the system:

- 1. Start the instrument (page 22).
- 2. Start the computer (page 24).
- 3. Check maintenance notifications in the Dashboard (page 28).
- 4. Check consumable status in the Dashboard (page 29).
- 5. Replenish consumables (page 31).

Set up and run: 1. Prepare the instrument (page 42). 2. Preheat the oven (page 42). 3. Check instrument status (page 53). 1. Create or import a plate (page 43). 2. Assign plate contents (page 46). Quick Start a run (page 55). 3. Print the plate layout (page 50). 4. Prepare and load sample plates (page 51). 1. Load plates for run and create the injection list (page 56). 2. Review and modify the injection list in Preview Run (page 59). 3. Start the run (page 61). 4. Monitor the run (page 61), check sequence or sample quality and specify re-injections (page 62). Review fragment/HID results: Review sequencing results: 1. Review sequence quality (page 81). 1. Review sample quality (page 89). 2. Specify re-injections (page 95). 2. Specify re-injections (page 85). 3. Review quality reports (page 85). 3. Review quality reports (page 95). 4. Export sequencing results (page 87). 4. Export sizing results (page 96). (Optional) print or save (.pdf) calibration and performance check reports to save with results: Spatial calibration (page 99) • Spectral calibration (page 103). .

- Sequencing install standard performance check (page 122).
- Fragment or HID install standard performance check (page 122).
- Fragment of Fild Install standard performance check (page 15

# Start the instrument

1. Verify that the instrument is connected to the appropriate power supply.

**CAUTION!** Do not unpack or plug in any components until the Applied Biosystems service representative has configured the system for the proper operating voltage.

See the *Applied Biosystems 3500 Series Genetic Analyzer Site Preparation Guide* (4401689) for details.

**Note:** The purpose of the Site Prep Guide is to help you prepare your site for installation of the 3500 or 3500xL analyzer. For specific details about your system, please refer to this user guide.

**IMPORTANT!** Do not rename the computer after the 3500 Series Data Collection Software has been installed. The instrument computer has been assigned a unique name. Changing the name may cause the 3500 Series Data Collection Software to malfunction.

- 2. Inspect instrument interior. Ensure that:
  - a. The oven door is closed.
  - b. No objects are left inside the instrument.

**IMPORTANT!** Misplaced objects left inside the instrument can cause damage.

- 3. Close instrument door.
- 4. Turn on the instrument. Press the power on/off button on the front of the instrument and wait for the green status light to turn on.



a. Press the Tray button on the outside of the instrument to bring the autosampler to forward position. Wait until the autosampler stops at the forward position.

**Note:** When the door is open, the yellow status light blinks while the instrument performs self-check and the autosampler adjusts.

b. Check the instrument status. Ensure the green status light is on and not flashing before proceeding. The table below explains the status indicator lights for the instrument.

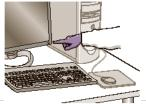
Indicator	Status		
All lights off	Instrument off		
Green light	Operational (awaiting run)		
	Pause run, terminate run, stop injection button (in SW) pressed by user.		
	<b>Note:</b> You can only abort an injection when the green light is flashing, not when it is solid green.		
Green light (blinking)	Operational (Run in progress)		
Amber light (blinking)	Power-up self-test in progress		
	Run paused		
	Door open		
	Run failure that doesn't require restart of instrument		
Amber light	Standby		
Red light	Self-test failed		
	Instrument failure		
	Requires a restart of the instrument and computer		

# Start the computer

1. Power on the computer.

- 2. Power on the monitor.
- 3. In the Log On to Windows dialog box:
  - a. Enter the user name.
  - b. If applicable, enter a password.





**Note:** If the computer is connected to a network, you do not need to log on to the network before starting the instrument.

c. Click OK. Wait until the computer finishes booting.

**IMPORTANT!** The status icon, on the right lower-corner of your screen, shows when the 3500 Server Monitor is active by displaying the icon shown here.



**IMPORTANT!** Do not close this icon. Doing so will prevent proper functioning of the software.

### Log on to Windows

Follow the prompts to log on to the Windows operating system.

## Launch the application

Step one: Launch<br/>the DaemonIf the Daemon does not start automatically, launch the DaemonStart ▶ Programs ▶ Applied Biosystems ▶ 3500 ▶ Daemon

👅 Daemon

Note: It will take approximately 15 seconds for Daemon to populate.

Step two: Launch the Server Monitor	If the Server Monitor does not start automatically, launch the Server Monitor: <b>Start &gt; Programs &gt; Applied Biosystems &gt; 3500 &gt; Server Monitor</b> Server Monitor	
	Note: It will take approximately 2 minutes for the Server Monitor to set up. During this time, you will see the status icon transition from a red circle, with an X in the middle (indicating that not all 3500 services are loaded) to the shape of an hour-glass on your desktop, next to the clock.	
	When Server Monitor set up is complete, the icon in the shape of an hour-glass will disappear and a checkmark icon appears indicating that the 3500 Server Monitor has started and all 3500 services loaded.	
Step three: Launch the 3500 application	Launch the application: Start → Programs → Applied Biosystems → 3500 3500	
Splash screen	After you launched the 3500 application, the 3500 Series Data Collection Software splach screen appears. This screen will remain active for a few seconds until the 3500	

splash screen appears. This screen will remain active for a few seconds until the 3500 Log In dialog box opens.



After the 3500 Series Data Collection Software splash screen disappears, one of the following occurs:

- ٠ The Dashboard is displayed (go to "Check system status in the Dashboard" on page 26)
- The Login dialog box is displayed (go to "Log In" on page 26) ٠

### Log In

Security, Audit, The Security, Audit, E-Signature (SAE) module is an optional component of the 3500 Series Data Collection Software. Researchers have the option to purchase this feature and enable/disable the functionality for SAE. If the SAE feature is enabled, see Chapter 7, Use Security, Audit, and E-Sig Functions (SAE Module) for user configurations.

After the 3500 Series Data Collection Software splash screen disappears, log in from the Dashboard:

1. Enter the User Name and Password in the 3500 Log In dialog box.

🐸 3500 Log In	$\mathbf{X}$
<b>3500 Log In</b> Provide your user name and password to login.	
User Name: Password:	
Reset	OK Cancel

2. Click OK.

The 3500 Series Data Collection Software splash screen re-appears. This screen will remain active for a few seconds and the 3500 Series Data Collection Software opens.

The 3500 Series Data Collection Software launches and the Dashboard appears.

**IMPORTANT!** If you accidentally close any of the services (via 3500 Server Monitor), the system will not work. To open a closed service, place the cursor on the status icon, click the right-mouse button, go to Services, and click the service that is closed.

## Check system status in the Dashboard

### Dashboard, a quick glance

The first screen that is displayed when you start the 3500 Series Data Collection Software is the Dashboard (Figure 4).

The Dashboard displays gauges, instrument information, consumable information, and maintenance notifications that provide a quick overview of the usage of each consumable and the status of the instrument.

Consumable containers include radio frequency identification (RFID) tags that identify the consumable and allow the software to monitor the number of runs or days remaining, the number of days on the instrument, the expiration date, lot number and part numbers.

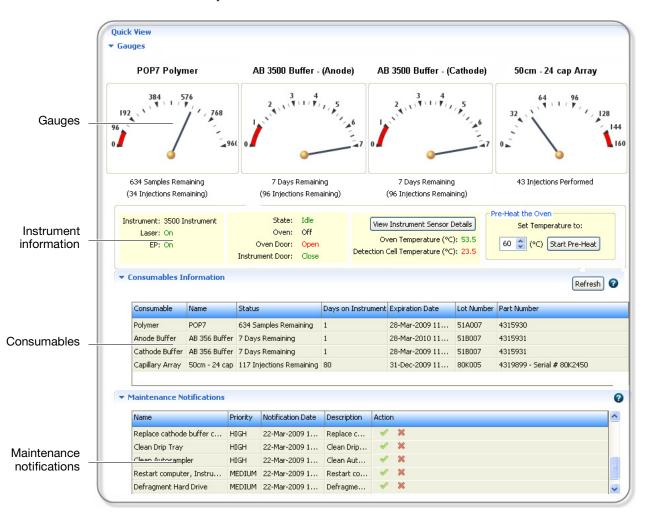


Figure 4 Dashboard

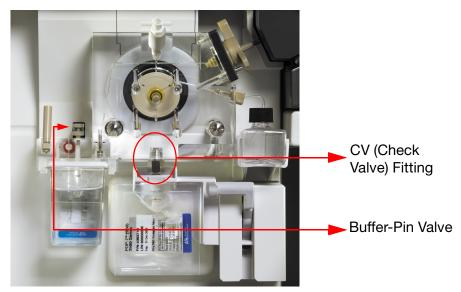
### Check maintenance notifications

The Maintenance Notification section displays reminders for the tasks scheduled in the maintenance calendar (see "Use the maintenance calendar" on page 232). You can set the time to trigger maintenance notifications in Preferences (see "Set general preferences" on page 33).

1. Review the Maintenance Notifications pane.

Name	Priority	Notification Date	Description	Action	1
Perform Performance Check	HIGH	28-Jan-2009 12:00:00 AM	Performance Check	<ul> <li>Image: A second s</li></ul>	ж
Clean Drip Tray	HIGH	28-Jan-2009 12:00:00 AM	Clean Drip Tray	-	ж
Clean Autosampler	HIGH	28-Jan-2009 12:00:00 AM	Clean Autosampler	×	х
Replace Reservoir Septa	HIGH	28-Jan-2009 12:00:00 AM	Replace Reservoir Septa	×	ж
Wash Pump Trap	HIGH	28-Jan-2009 12:00:00 AM	Wash Pump Trap	1	х

- 2. Perform any scheduled maintenance tasks, then click in to mark it as complete, (or click in to mark it as dismissed if you do not perform the task). Actions are recorded in the Notifications log (for more information, see "Review the Maintenance Notifications Log" on page 257).
- 3. Perform any daily, monthly, or quarterly maintenance tasks that are not listed in the Maintenance Notifications pane (see Chapter 8, Maintain the Instrument).
- 4. Inspect the instrument interior. See "Start the instrument" on page 22.
  - a. If you see any spills, clean immediately.
  - b. If you see any leaks and dried residue around the Buffer-Pin Valve, check valve, and array locking lever. If leaks persist, contact Applied Biosystems.



### Check consumable status

**IMPORTANT!** The Days Remaining for buffers updates only when you click Refresh or start a run. As part of daily startup, click **Refresh** to update consumable status.

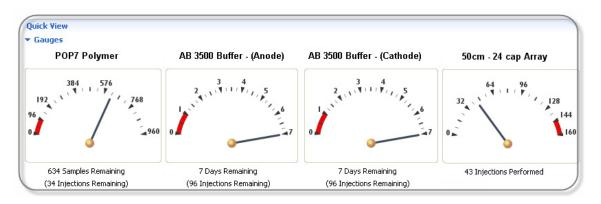
1. Click **Refresh** to update consumable status.

The Consumables pane displays expiration dates and lot numbers (read from the RFID tags on the consumable containers).

Consumables Information     Refresh						Refresh 🕜
Consumable	Name	Status	Days on Instrument	Expiration Date	Lot Number	Part Number
Polymer	POP7	634 Samples Remaining	1	28-Mar-2009 11	51A007	4315930
Anode Buffer	AB 356 Buffer	5 Days Remaining	1	28-Mar-2010 11	51B007	4315931
Cathode Buffer	AB 356 Buffer	5 Days Remaining	1	28-Mar-2010 11	51B007	4315931
Capillary Array	50cm - 24 cap	117 Injections Remaining	80	31-Dec-2009 11	80K005	4319899 - Serial # 80K2450

2. Check the consumables gauges for the number of injections, samples, or days remaining for a consumable. Table below lists specifications for each consumable.

When <10% of the specified use of the consumable remains, the gauge moves into the red warning range. The consumable also displays in red in the Consumables pane.



**IMPORTANT!** Applied Biosystems recommends that you add a maintenance notification to your calendar for polymer and buffer replacement. Set the notification to display two days before the polymer should be replaced.

Consumable		mable	On-instrument limits (the first limit met applies)	Notes	
Polymer <sup>‡§</sup>	8-cap	960 sample pouch	960 samples or 120 injections	Use within 7 days of installation on	
		384 sample pouch	384 samples or 60 injections	instrument.	
	24-cap	960 sample pouch	960 samples or 50 injections	<ul> <li>The software allows you to continue running past 7 days. However,</li> </ul>	
		384 sample pouch	384 samples or 20 injections	Applied Biosystems has verified the polymers for up to 7 days only on the instrument.	
Buffers	Buffers 8-cap		7 days or 120 injections	To ensure optimal buffer	
24-cap			7 days or 50 injections	performance, the software requires buffer replacement after 7 days.	
Capillary Array			160 injections	The software allows you to continue running after 160 injections. However, Applied Biosystems has verified the arrays for up to 160 injections.	

The Polymer Sample Counter decrements only for wells that contain sample, but the Polymer Injection counter decrements for each injection, regardless of whether all wells contain sample. The sample limit and the corresponding injection limit may not coincide. Note that the initial injection limit is higher than the initial sample limit.

Example: 960 sample pouch on 24-cap:

If all wells contain sample for all injections: 960/24 = 40 injections.

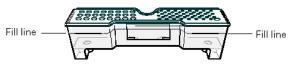
If all wells do not contain sample for all injections: 960/<24 = 40+ injections, up to a maximum of 50 injections and a maximum of 960 samples.

A polymer pouch includes additional volume to accommodate the volume used during installation and by wizards. However, excessive use of wizards reduces the number of remaining samples and injections, based on how many times specific wizards are run. For example, if you run the total bubble remove option in the Remove Bubbles wizard more than four times or run other wizards excessively, including multiple pouch installations, the number of remaining samples and injections is reduced.

including multiple pouch installations, the number of remaining samples and injections is reduced. § Ambient temperature must be in the range of 15 °C to 25 °C POP-6<sup>™</sup>. Sustained use at higher temperatures may result in shorter read lengths than specified.

## Check buffer fill levels

Check the fill levels on buffers. Verify that buffer level is at the top of the fill line and check that seal is intact.



**IMPORTANT!** Do not use if the buffer level is too low or the seal has been compromised. Ensure that the buffer level is at or above the fill line and the seals is intact.

## **Replenish consumables**

As needed, see:

- "Replenish polymer" on page 245.
- "Change polymer type" on page 247.

**IMPORTANT!** Wear gloves while handling polymer, the capillary array, septa, or CBC.

- "Change the anode buffer container (ABC)" on page 237.
- "Change the cathode buffer container (CBC)" on page 238.
- "Fill capillary array with fresh polymer" on page 251
- "To change the capillary array" on page 252.

Go to Chapter 3, "Set Up and Run" on page 41.

# Set preferences

## **Overview**

Preferences are user-definable default settings. To access the Preferences dialog box, select Preferences in the toolbar. You can optionally set any or all preferences.



Je Preferences	
type filter text not used	System 🗘 🗢 🔿 🗠
System     Date Format     Date Format     Date Format     Scheduler Preference     Sequencing Settings     Deter     Spectral Calibration     Deter     Plate Setup     Reports Settings     Run Setup     Sequencing Settings	Select one of the sub-categories to set its properties. You may choose to reset all system preferences to the factory installed defaults. If your user permissions enable you to do so, you may import or export the system preferences.
<ul> <li>Sequencing Settings</li> <li>Trace</li> <li>Trace Print</li> <li>Trace Quality</li> <li>Trace Quality Reports</li> </ul>	Import Export Restore Defaults Apply
	OK Cancel

Note: The "type filter text" field at the top of the dialog box is not used.

## System preferences

These settings apply to all users:

- Date format
- Instrument settings (instrument name)
- Scheduler preference (trigger time for maintenance notifications)
- Sequencing export settings
- Spectral calibration (number of allowed borrowing events)

### **User preferences**

These settings apply to all users if your system does not include the SAE module, but are saved individually per user if your system includes the SAE module:

**Note:** For information on the SAE module, see Chapter 7, "Use Security, Audit, and E-Sig Functions (SAE Module)" on page 197.

- Plate setup
- Reports settings
- Run setup
- Sequencing (review and report settings)

### Set general preferences

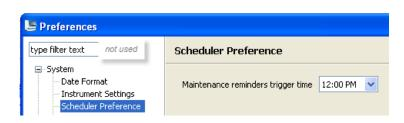
- **System** 1. In the Preferences dialog box, click the following items:
- preferences
- **Date Format** to set the date and time format for the software.



• **Instrument Settings** to set the instrument name (appears in the Dashboard, reports, file name conventions, instrument sensor details, view sequencing results).

Je Preferences	
type filter text not used	Instrument Settings
- System - Date Format - Instrument Settings	Instrument name 3500 Instrument

• Scheduler Preference to set the time to trigger maintenance notifications displayed in the Dashboard (see "Check maintenance notifications" on page 28).



• **Spectral Calibration** to decrease the number of allowed borrowing events for spectral calibration (see "What you see during a spectral calibration" on page 112).

🐸 Preferences	
type filter text not used	Spectral Calibration $\diamond \star \Rightarrow \star$
<ul> <li>System</li> <li>Date Format</li> <li>Instrument Settings</li> <li>Scheduler Preference</li> <li>Sequencing Settings</li> <li>Export</li> <li>Spectral Calibration</li> </ul>	Choose the maximum number of borrowing events for a 24-Cap array. 3

- 2. Click **Apply** to save the system preferences (see "System preferences" on page 32).
- **User preferences** 1. In the Preferences dialog box, click the following items as needed:
  - Plate setup to set the default settings for:
    - Plate type and attributes when you create a plate.
    - Plate type in the Open Plate dialog box.

블 Preferences	
type filter text not used	Plate Setup 🗢 🔹 🔶 🕤
⊕- System ⊡- User	Choose the default application type, polymer type, and capillary length to use when the instrument settings cannot be detected.
···· <mark>Plate Setup</mark> ···· Reports Settings ···· Run Setup	Application Sequencing V Polymer Type POP7 V Capillary Length 50 V
<ul> <li>Sequencing Settings</li> <li>Trace</li> <li>Trace Print</li> </ul>	Choose the default Assign Plate Contents View.
Trace Quality Trace Quality Reports	Assign Plate Contents View Plate 💌

**Reports settings** to set the default font and size reports.

Note: You can override this setting in each report view.

블 Preferences	
type filter text not used ⊪-System	Reports Settings These settings affect all reports
User Plate Setup Reports Settings Run Setup Sequencing Settings Trace	Font settings Select the font to be used in reports.
Trace Print Trace Quality Trace Quality Reports	

• **Run Setup** to set the default storage location for data files in file name conventions and results groups.

**Note:** You can override this setting in file name conventions and results groups.

Je Preferences		
type filter text not used	Run Setup	\$• € *
System     System     Vser     Plate Setup	Choose the default file location where sample files w D:\Applied Biosystems\3500\Data	Ill be stored. Browse
Run Setup Sequencing Settings		

2. Click Apply to save the user preferences (see "User preferences" on page 33).

Table and plot settings user preferences Users can also save user preferences while viewing tables and plots:

- **Table settings dialog box** Determines the columns displayed in a table and the order of the columns.
- Plot settings dialog box Determines the settings applied to plots.

### Set sequencing preferences

**Export (system preference)** Export preferences set the defaults for the file types to automatically export during a sequencing run. Exported files are stored in the same directory as the .ab1 files.

1. In the Preferences dialog box, click **Export** under System Sequencing settings to display the Export pane.

🔚 Preferences			
type filter text not used	Export		
	Select the file types to export to export the entire sequence File Type: *.annotation.txt *.phd.1 *.scf		check each file type and specify :k Apply.
Plate Setup Reports Settings		Entire Sequence	Post-trim Sequence Only
- Repetered Settings - Sequencing Settings - Trace - Trace Print	*.fsta	0	⊙
Trace Quality	- *,qual	0	۲
	*.seq	0	۲

2. Select the file types to export. Exported files are stored in the same directory as the .ab1 files.

File type	Description
*.annotation.txt	Information from the Annotation tab in the sequencing trace view such as data collection time, run time start finish
*.phd.1, *.scf	Sequencing files
*.fsta, *.qual, *.seq	Reference files – specify Entire Sequence or Post-trim Sequence Only

3. Click **Apply** to save the system preferences (see "System preferences" on page 33).

**Trace (user preference)** The Trace preference settings determine the default settings for color representation of nucleotide and quality value bars in the Trace View in View Sequencing Results.

1. In the Preferences dialog box, click **Trace** under User Sequencing settings to display the Trace pane.

Je Preferences		- 🗆 🗙
type filter text not used	Trace	י לף י
System     System     User     Plate Setup     Reports Settings     Run Setup     Sequencing Settings     Trace     Trace Print     Trace Quality	These settings affect the Trace panel.          NT Base Color       Pure Base QV Colors         NT Base Foreground Background       0         A       Image: Color Colors         G       Image: Color Colors	100
Trace Quality Reports	T Mixed Base QV Colors 0 10 15	100
	Mixed Preview	
	Level Market and the first state of the stat	
	CAWTACGAAKAAAGTCTGAAACTGTTG 10 20	AGAG 30

2. Specify the following settings.:

Setting	Description
NT (nucleotide) Base Color	Click an NT or Mixed base Foreground or Background color block, then select a color for the letter annotation or the highlight color for the letter annotation.
Pure Base and Mixed Base QV Colors	Sets the colors and ranges for pure and mixed base quality value indicators (QVs) displayed in the Trace View (the default settings are recommended):
	a. Click a pure base or mixed base color bar to select a new color.
	b. Place the mouse pointer over a slider, then drag to set a new range.
	Applied Biosystems recommends that you set the following ranges for QVs:
	<ul> <li>Pure bases: Low QV ≤ 15, Medium QV = 15 to 19, High QV = 20+ (default)</li> </ul>
	• Mixed bases: Low $QV \le 5$ , Medium $QV = 5$ to 10, High $QV > 10$ (investigate to determine the best range for your application)
	Note: The predicted probability of error for a basecall is high QV >10.

3. Click **Apply** to save the user preferences (see "User preferences" on page 34).

### Trace Print (user preference)

Trace Print preferences determine settings for sequencing trace reports.

1. In the Preferences dialog box, click Trace Print under User Sequencing settings to display the Trace Print pane.

🔄 Preferences	
type filter text not used	Trace Print
- System - User - Plate Setup - Reports Settings - Run Setup - Sequencing Settings - Trace - Trace Print - Trace Quality - Trace Quality Reports - Tr	These settings affect the printed Trace Reports Select data to be printed for each trace Analyzed data  Entire sequence Post-trim sequence only Annotation Sequence Select Settings Panels per page  Points per panel 1000 Set Y-Scale for analyzed data  Individual Uniform with a maximum sc 4000

- 2. Specify the type of trace data, specific print settings, and Y-Scale preference to display in the Trace Report.
- 3. Click Apply to save the user preferences (see "User preferences" on page 34).

**Trace Quality** Trace Quality preferences control the quality ranges for: (user preference)

- **QC report** Trace Score and CRL
- **Plate report** Trace Score
- 1. In the Preferences dialog box, click Trace Quality under User Sequencing settings to display the Trace Quality pane.

type filter text not used	Trace Quality $\diamond \star \Rightarrow$
System User Very Reports Settings Run Setup Sequencing Settings Trace Trace Print Trace Quality Trace Quality Reports	These settings define thresholds for the Trace Quality Reports. Trace scores and contiguous read length settings can also contribute colors used within the Trace Quality Reports. To change these settings, drag the sliding bar to the desired setting. Alternatively, right click on the bar and enter the desired setting. Trace Score Color and Threshold Setting The Trace Score is the average of basecall quality values for bases in the clear range. 0 15 30 100 Contiguous Read Length Color and Threshold Setting Full Sequence Drop In Quality Yalues Contiguous Read Length (CRL) The Contiguous Read Length (CRL) is the longest uninterrupted stretch of bases with quality higher than a specified limit. It is the length of the sequence that remains after trimming only using quality values. Evaluation of the quality of each base is determined using the quality value of adjacent bases within a specified window. 0 10 300 1000

- a. Click a color bar to select a new color.
- b. Place the mouse pointer over a slider, then drag to set a new range.



3. Click **Apply** to save the user preferences (see "User preferences" on page 34).

### Trace Quality Report (user preference)

Trace quality Report preferences determine the content and formatting used in QC, Plate, Trace Score, CRL, QV20+, and Signal Strength reports.

1. In the Preferences dialog box, click **Trace Quality Report** under User Sequencing settings to display the Trace Quality Report pane.

🐸 Preferences	
type filter text not used	Trace Quality Reports $\Leftrightarrow \neg \Rightarrow \neg$
<ul> <li>System</li> <li>Date Format</li> <li>Instrument Settings</li> <li>Scheduler Preference</li> <li>Sequencing Settings</li> <li>Export</li> <li>Spectral Calibration</li> <li>User</li> <li>Plate Setup</li> <li>Reports Settings</li> <li>Run Setup</li> <li>Sequencing Settings</li> <li>Trace</li> <li>Trace Print</li> <li>Trace Quality</li> <li>Trace Quality Reports</li> </ul>	These settings affect all reports         Sort data based on         This setting applies to the Trace Score Report, CRL Report, QV20+ Report, and Signal Strength Report.            • Run Name         • Capillary Number          Signal based on         This setting applies to the QC Report, and Signal Strength Report.            • Average Raw Signal Intensity         • Average Raw Signal to Noise Ratio          Display well image by            • Wider thumbnail without the file name         • Smaller thumbnail with the file name         • Smaller th

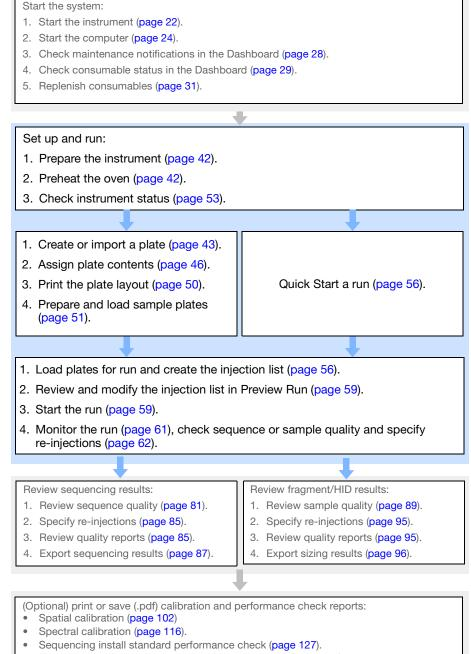
2. Specify the following settings.:

Setting	Description				
Sort data	Sort data in Trace Score, CRL, QV20+, and Signal Strength reports based on:				
	Run Name				
	Capillary Number				
Signal based on	Base signal in QC and Signal Strength reports based on:				
	Average Raw Signal Intensity				
	Average Raw Signal to Noise Ratio				
Display well	Specify the thumbnail option for Plate reports:				
image by	Wider thumbnail without file name				
	Smaller thumbnail without file name				

3. Click Apply to save the user preferences (see "User preferences" on page 34).

# Set Up and Run

## Workflow



# Prepare the instrument

- 1. In the Dashboard, check consumable status (page 29). Ensure that:
  - Consumables are not expired
  - Adequate buffer levels are at the fill lines.
- 2. Set the oven temperature, then click **Start Pre-heat**:
  - 60 °C POP-7<sup>TM</sup> and POP-4<sup>TM</sup> polymers
  - 50 °C POP-6<sup>TM</sup> polymer

Pre-heat the oven and detection cell while you prepare for a run (detection cell temperature is set by the software). Preheating helps mitigate subtle first-run migration rate effects. The preheat function automatically turns off after 2 hours of instrument inactivity.

Applied Biosystems recommends that you pre-heat the oven for at least 30 minutes before you start a run if the instrument is cold.

3. Check the pump assembly for bubbles and run the Remove Bubble wizard if needed (see page 251).

## Create a plate

**Note:** If you are running a stand-alone version of the 3500 Series Data Collection Software (a version that is not installed on the instrument computer), you can create plates, then export them for use on the instrument computer.

# Create a plate from a template

The software includes factory-provided plate templates that you can use as a starting point to create a plate (you can also create your own plate templates). In addition to pre-defined plate parameters, a plate template can also contain a list of the appropriate assays, file name conventions, and results groups for an application. For more information, see "Create a plate template" on page 75.

 In the Dashboard, click Create Plate From Template to display the Open Plate Template from Library dialog box.



<b>U </b> 0 p	en Plate Template From	Library	×					
	Instructions Select row from table and click on "Open" button.							
Fil	ter: Sequencing	Sea	rch: All So Clear 😮					
	Plate Name	Туре	Description					
1	AR BDx_Rapid_Seq	Sequencing	For use with samples purified with BigDye XTermin					
2	🛝 📝 Std_Seq-POP7	Sequencing	For read lengths of 850 bp or greater and a run til					
3	🛝 📝 Short_Read_Seq	Sequencing	For read lengths of 300 bp and a run time of 30 m					
4	AR Std_Seq_xL-POP6	Sequencing	For read lenghts of 600 bp or greater and a run ti					
5	AR Std_Seq_xL-POP7	Sequencing	For read lengths of 850 bp or greater and a run ti					
6	ARS 🔀 Fast_Seq_xL-POP7	Sequencing	For read lengths of 700 bp and a run time of 65 m					
7	ARB BDx_Fast_Seq_xL	Sequencing	For use with samples purified with BigDye XTermin					
8	ARS BDx_Std_Seq-POP7	Sequencing	For use with samples purified with BigDye XTermin					
<	AS POR CHI Can DODA	Sequencing	Sor use with complex purified with RieDue VTermin					
	īlose		Open					

- 2. (Optional) Filter the templates listed:
  - a. Select a template type.

Filter:	Sequencing	~	
	All		
	Sequencing		
Pla	Fragment Mixed (Seq + Fi HID	rag)	
Search:	30	All	GO Clear
		Plate Name	
		Туре	
Desc	ription	Barcode	~

**Note:** You can set the default plate type for this filter in Preferences. See "Specify the default plate type for the Open Plate dialog box" on page 76.

- b. Find templates by selecting an attribute, entering the text to search for, then clicking **Go**. (Click **Clear** to clear the field and enter different search criteria).
- 3. Select the template, then click **Open**.
- 4. In the Define Plate Properties screen, select the plate type.

📖 New Plate 🔻 🔛 Open Plate 🔻 🔛 Save Plate 🔻 🛄 Close Plate
Plate Details
* Name: Enter plate name
* Number of Wells: 💿 96 🔵 96-FastTube 🔵 384
* Plate Type: Sequencing
* Capillary Length: 50 🗸 cm
* Polymer: POP7 🗸

- **96** Supports 96-well standard reaction plate. 8-strip standard tubes are also supported with appropriate retainers.
- **96-Fast Tube** Supports 96-well Fast reaction plate. 8-strip fast tubes are also supported with appropriate retainers.
- 5. Set remaining plate properties, then select Save.
- 6. Click Assign Plate Contents, then go to "Assign plate contents" on page 46.

**Import a plate** 1. Do either of the following:

- Create a plate on another 3500 Series Data Collection Software system, then export (see "Import and export a plate" on page 75).
- Create a plate import file (see "Create a plate import file" on page 74).

- 2. Access the Assign Plate Contents screen: Click the Main workflow arrow ( , in the Dashboard, then select Assign Plate Contents in the navigation pane.
- 3. Click **import**, then select the plate import file.
- 4. Click Assign Plate Contents.



# Assign plate contents

You assign the following information to the wells in a plate before you can run the plate:

- Sample names and sample types (required) Identifies the well positions of each sample for data collection and processing.
- Assay (required) Specifies the parameters that control data collection and primary analysis (basecalling or sizing). All named wells on a plate must have an assigned assay. For more information on assays, see "Assays library" on page 147.
- Filename convention (optional) Specifies file naming. For more information on assays, see "File name conventions library" on page 151.
- **Results group** (optional) Specifies sample data file storage. For more information on assays, see "Result group library" on page 155.

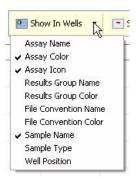
Before you assign plate contents

- 1. Access the Assign Plate Contents screen (Figure 5 on page 47) from:
  - The Define Plate Properties screen by clicking **Assign Plate Contents** (described above).



- The navigation pane by selecting **Assign Plate Contents** in the navigation pane.
- The Dashboard by clicking the Main workflow arrow ( , then selecting Assign Plate Contents in the navigation pane.
- 2. Create a plate. Select one of the following topics:
  - "Create a new plate" on page 144
  - "Create a plate from a template" on page 43
  - "Import a plate" on page 44
  - Or select Open Plate > Edit Existing Plate
- 3. Click **Show In Wells** to specify the attributes to display in wells.

Figure 5 on page 47 shows the Plate View of the Assign Plate Contents screen.

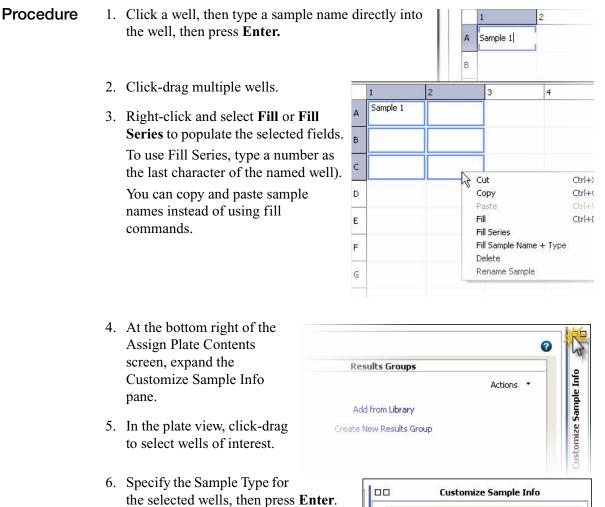


											— Show well attributes
Show In Well	; • Sele	ct Wells 🔻	Array	/ Selection	Row	E Colun	nn 🛛 🔝 2	loom In 🚦	🔄 Zoom (	Dut 🔣	
1 2	3	4	5	6	7	8	9	10	11	12	
A											
B											Name samples
C D											
E											
F											
G											
н											
				-							Assign assays, file
Name:				Ba	arcode:						<ul> <li>name conventions, and results groups</li> </ul>
Name:					arcode:	Cus	tomize S	ample Inf	0		
Name:	Assa				D D Property	,	tomize S	· ·	0		and results groups
Vame:	Assa	γs	Actions	0	Property I 1: Sa	/ mple ample Nam	Value e		o		and results groups
	Assa 1_Assay_xL-POF		Actions	0	Property I: Sa S 2: Cu	/ mple ample Nam ample Type istom	Value e Sampl		o 		and results groups
				0	Property I: Sa S S 2: Cu L	/ mple ample Nam ample Type	Value e Sampl d Fie		0		Assign sample types and

Figure 5 Assign Plate View of the Assign Plate Contents screen

## Name samples and assign sample types in the plate view

This section provides one way to name samples and assign sample types. For other ways to name samples, see "Use the Plate View" on page 70 and "Use the Table View" on page 71.



 (Optional) Specify User Defined Fields and Comments. User Defined Fields contain additional attributes you can assign to a plate and are displayed only in Table View.

	Custor	nize Sample Info	
Pr	operty	Value	
	1: Sample		
	Sample Name		
	Sample Type	Sample	V
	2: Custom	Sample	^
	User Defined	Positive Control	
	User Defined	Allelic Ladder	
	User Defined	HiDi	~
	User Defined		
	User Defined		
-	3: Misc		
	Comments		

- 8. (Optional) For sequencing assays, specify amplicon and specimen.
- 9. Repeat to assign the Sample Type for all named wells.
- 10. Go to "Assign assay, file name convention, and results group in the Plate View" on page 49.

Property	Value
🖃 1: Sample	
Sample Name	
Sample Type	Sample
2: Custom	
User Defined Field 1	
User Defined Field 2	
User Defined Field 3	
User Defined Field 4	
User Defined Field 5	
🗐 3: Analysis	
Amplicon	
Specimen /	
E 4: Misc	

**Note:** For HID applications, include the well position in the allelic ladder sample names. Well position is needed to identify the position of allelic ladder samples during re-injection.

## Assign assay, file name convention, and results group in the Plate View

**Note:** If an assay, file name convention, or results group is not listed for the plate, go to "Add assays, file name conventions, and results groups to a plate" on page 73.

- 1. Select the wells for which to specify an assay.
- 2. Enable the checkbox next to the assay name to assign it to the selected wells.

Assays			
		Actions	•
🕑 🛃 BDx Fast Seq Assay xL - P	2	X	

**Note:** To normalize fragment analysis or HID data, select an

assay that contains a sizecalling protocol or a QC protocol that specifies a normalization size standard.

3. (Optional) Repeat for file name conventions and results group.

File Nam	e Conventions	Results Groups	
	Actions 🔻		Action
O My FNC	2	💿 🧹 My Sequencing Results Gro 📝	X
12		12	

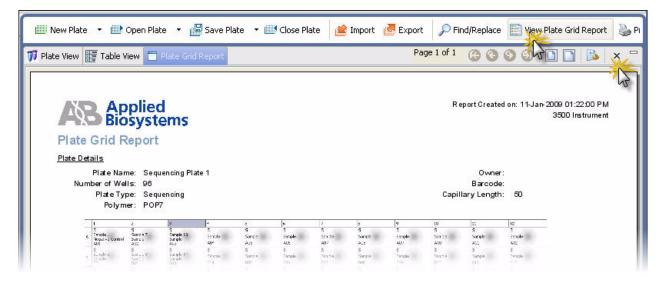
- 4. Select Save Plate.
- 5. Go to "Print the plate layout" on page 50.

How file location<br/>in file name<br/>conventions and<br/>results groups<br/>workIf you do not specify a file name convention, data files are named in this format:<br/><sample name>\_<well>.If you do not specify a results group, files are stored in the location specified in the<br/>file name convention or in Preferences > User > Run (see "User preferences" on<br/>page 34).

If you specify both a file name convention and a results group, files are stored in the location specified in the results group.

## Print the plate layout

1. In the Assign Plates for Run screen, click View Plate Grid Report.



**Note:** A 384-well report displays the plate layout in four quadrants on four pages.

- 2. Select Print Preview or Print as needed.
- 3. To save the report electronically (.pdf), print the report and select **CutePDF** Writer as the printer.
- 4. Close the report.



5. Go to "Prepare and load sample plates" on page 51.

# Prepare and load sample plates

**IMPORTANT!** Do not use warped or damaged plates.

# Capillary-to-plate mapping

The capillary-to-plate mapping for the default injection order is shown below. If you change the injection order in the injection list, mapping differs from the examples shown below.

- **96** Supports 96-well standard reaction plate. 8-strip standard tubes are also supported with appropriate retainers.
- **96-Fast Tube** Supports 96-well Fast reaction plate. 8-strip fast tubes are also supported with appropriate retainers.

#### 8-capillary: 96-well plate

	Cap		3	4	5	6	7	8	9	10	11	12
Α	1	1	3	4	5	6	7	8	9	10	11	12
В	2	- 1	3	4	5	6	7	8	9	10	11	12
С	3.	- 1	3	4	5	6	7	8	9	10	11	12
D	1	2	3	4	5	6	7	8	9	10	11	12
E	1	2	3	4	5	6	7	8	9	10	11	12
F	1	2	3	4	5	6	7	8	9	10	11	12
G	1	2	3	4	5	6	7	8	9	10	11	12
н	1	2	3	4	5	6	7	8	9	10	11	12

#### 24-capillary: 96-well plate

	1	2	3	4	5	6	7	8	9	10	11	12
Ca	p 1	2	3	2	2	2	3	3	3	4	4	4
	4	5	6	2	2	2	3	3	3	4	4	4
C	1	1	1	2	2	2	3	3	3	4	4	4
D	1	1	1	2	2	2	3	3	3	4	4	4
Е	1	1	1	2	2	2	3	3	3	4	4	4
F	1	1	1	2	2	2	3	3	3	4	4	4
G	1	1	1	2	2	2	3	3	3	4	4	4
н	1	1	1	2	2	2	3	3	3	4	4	4

#### 8-capillary: 384-well plate

Not supported on the 3500 Dx Genetic Analyzers (8-capillary)

Ca	D				
-	1	2	3	4	5
A	1	3	2	3	3
В	2	4	2	4	2
C	4	3	5	3	6
D	2	4	2	4	2

#### 24-capillary: 384-well plate

	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24
A	1	3	1	3	1	3	5	7	5	7	5	7	9	11	9	11	9	11	13	15	13	15	13	1
в	2	4	2	4	2	4	6	8	6	8	6	8	10	12	10	12	10	12	14	16	14	16	14	10
С	1	3	1	3	1	3	5	7	5	7	5	7	9	11	9	11	9	11	13	15	13	15	13	1
D	2	4	2	4	2	4	6	8	6	8	6	8	10	12	10	12	10	12	14	16	14	16	14	1
E	1	3	1	3	1	3	5	7	5	7	5	7	9	11	9	11	9	11	13	15	13	15	13	1
F	2	4	2	4	2	4	6	8	6	8	6	8	10	12	10	12	10	12	14	16	14	16	14	1
G	1	3	1	3	1	3	5	7	5	7	5	7	9	11	9	11	9	11	13	15	13	15	13	1
н	2	4	2	4	2	4	6	8	6	8	6	8	10	12	10	12	10	12	14	16	14	16	14	1
I	1	3	1	3	1	3	5	7	5	7	5	7	9	11	9	11	9	11	13	15	13	15	13	1
J	2	4	2	4	2	4	6	8	6	8	6	8	10	12	10	12	10	12	14	16	14	16	14	1
К	1	3	1	3	1	3	5	7	5	7	5	7	9	11	9	11	9	11	13	15	13	15	13	1
L	2	4	2	4	2	4	6	8	6	8	6	8	10	12	10	12	10	12	14	16	14	16	14	1
М	1	3	1	3	1	3	5	7	5	7	5	7	9	11	9	11	9	11	13	15	13	15	13	1
N	2	4	2	4	2	4	6	8	6	8	6	8	10	12	10	12	10	12	14	16	14	16	14	1
0	1	3	1	3	1	3	5	7	5	7	5	7	9	11	9	11	9	11	13	15	13	15	13	1
P	2	4	2	4	2	4	6	8	6	8	6	8	10	12	10	12	10	12	14	16	14	16	14	1

### Allelic ladder run requirements

Applied Biosystems recommends that you inject one allelic ladder for each set of 24 samples:

- 8-capillary instruments One allelic ladder per 3 injections
- 24-capillary instruments One allelic ladder per 1 injections

Allelic ladders that are injected under the same conditions are recommended to accurately genotype samples in the secondary analysis software (GeneMapper<sup>®</sup> *ID-X* Software v1.2 or later).

**IMPORTANT!** Variation in laboratory temperature can cause changes in fragment migration speed that can, in turn, cause sizing variation. Applied Biosystems recommends the frequency of allelic ladder injections described above to account for normal variation in fragment migration speed. However, during internal HID validation studies, verify the required allelic ladder injection frequency to ensure accurate genotyping of all samples in your laboratory environment.

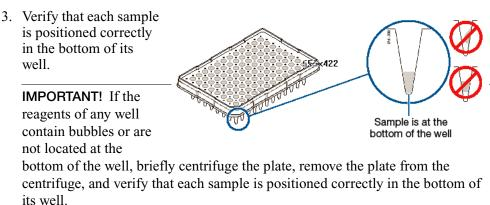
### Results group for one allelic ladder per run folder

For a 24-capillary instrument, create a results group that specifies an injection folder, then select this results group for all injections on the plate.

For an 8-capillary instrument, create one results group for each set of three injections on the plate (each results group specifies a results group name folder). For more information, see "Results group example 2: store one allelic ladder per run folder (8-capillary instruments)" on page 161.

### Prepare sample plates

- 1. Pipette samples into the plate according to the plate layout (see "Print the plate layout" on page 50).
- 2. Briefly centrifuge the plate.



4. Store the plate on ice until you prepare the plate assembly and load the plate in the instrument.

### Prepare the plate assembly

**IMPORTANT!** Prepare the plate assembly on a clean, level surface. Do not heat plates that are sealed with septa.

1. Align the holes in the septa strip with the wells of the plate, then firmly press downward onto the plate. 2. Place the sample plate into the plate base.

**IMPORTANT!** Make sure to use the correct plate base for standard plates versus 8-tube strips and fast plates. Using the wrong plate base may affect performance.

- Plate retainer Plate with septa strip Plate base
- 3. Snap the plate retainer (cover) onto the plate, septa, and plate base.
- 4. Verify that the holes of the plate retainer and the septa strip are aligned. If holes are not aligned, re-assemble and then assemble the plate assembly.

**IMPORTANT!** The array tips will be damaged if the plate retainer and septa strip holes do not align correctly.

## Load the plate in the instrument

1. Place the plate in the autosampler with the labels facing you (or the instrument door) and the notched corner of the plate in the notched corner of the autosampler.



2. Close the instrument door to re-initialize the instrument.

### Check instrument status

Check instrument status in the Dashboard. Temperatures are displayed in red as they warm to the set-points. When temperatures are at the set point they are displayed in green. Temperatures may fluctuate slightly when they reach the set point as they stabilize.

Applied Biosystems recommends that you pre-heat the oven for at least 30 minutes before you start a run if the instrument is cold. Pre-heating mitigates subtle first-run migration rate effects. (If you start the run when red indicators are shown, the run does not start until all indicators are green.)

Instrument: 3500 Instrument State: Idle Laser: On Oven: Off EP: On Oven Door: Open Instrument Door: Close	View Instrument Sensor Details Oven Temperature (°C): 53.5 Detection Cell Temperature (°C): 23.5	Pre-Heat the Oven Set Temperature to: 60 📚 (°C) Start Pre-Heat
--	--	--

## Link the plate

1. In the Assign Plates for Run screen, click Link Plate for Run.

📕 3500 Data C	Collection Softwar	е					
Dashboard	Edit Library Maint	tenance	Tools 🔹	Manage	<ul> <li>Prefer</li> </ul>	ences Hel	p 🔻 Logou
Plate Name:Seq		. 📖 Ne	ew Plate 🔹	🛄 Ор	en Plate	🕶 🔡 Sav	e Plate 🔻
AB Bid	oplied osystems	📆 Plate	e View 🕎	Table View	1		-
Setup Defi	ne Plate Properties	2	Show In V	/ells 🔻	- Sele	ct Wells 🔻	Arra
As	sign Plate Contents		1	2	3	4	5
🌄 Run Instrume	ent Load Plates for Run Preview Run Monitor Run	5 Nar	Sequencin		1		9 9; 9; 9; 9;
	ults Sequencing Results agment/HID Results				(2) Assi	Pr	:
	A COL			Link Pl	ate for Ru		*Mi

2. Go to "Load plates for run and create the injection list" on page 56.

Note: By default, the plate in position A is selected.

### Quick Start a run

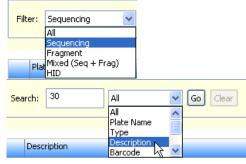
You can start a run in the Dashboard by selecting a plate with plate contents already assigned.

Load the plate in the instrument before proceeding (see "Load the plate in the instrument" on page 53).

1. In the Dashboard, click **Quick Start Run** to display the Select Plate from Library dialog box.

nstr				
	uctions			
Selec	t row from table and click o	n "Link Plate" butto	n.	
Filt	er: Sequencing	× 9	earch: All	
	Plate Name	Туре	Description	~
		Type	Description	
1	3-07-09 Run 1	Sequencing		
2	3-07-09 Run 2	Sequencing		~
<	ш			>
_	lose		Link	

- 2. (Optional) Filter the plates listed:
  - a. Select a plate type (you can set the default plate type in Preferences, see "Specify the default plate type for the Open Plate dialog box" on page 76).
  - b. Find a plate based on an attribute by selecting an attribute, entering the text to search for, then clicking Go. (Click Clear to clear the field and enter different search criteria).



»

Run

- 3. Select a plate, then click Load Plate.
- 4. Click Start Run from the Load Plates on Run Screen.

**IMPORTANT!** It takes, approximately, 10 seconds for the instrument to initialize after the instrument door is closed. Do not start a run until the instrument status light is green.

### Load plates for run and create the injection list

Load the plate in the instrument (see "Load the plate in the instrument" on page 53) and link the plate ("Link the plate" on page 54) before proceeding.

- 1. Access the Load Plates for Run screen (Figure 6 on page 56) from:
  - The Assign Plate Contents screen by clicking Link Plate for Run.



- The navigation pane by selecting Load Plates for Run in the navigation pane.
- The Dashboard by clicking the Main workflow arrow **(**, then selecting Load Plates for Run in the navigation pane.

	In Information	there has an head on the state							
		Name by entering text							
_		09-03-08-13-28-46-9	03	Connection St	atus: Connected		User Name: Administrat	or	
Pla	ates on Instrume	nt							
P	late A (96 wells)		Link Plate	Inlink Plate B			Link Plate Unlink	Recent Plates Re	ecent Runs
								Name	Date Modified
	Name: 3-09 Run	2						3-09 Run 2	08-Mar-200
	Type: HID			1				3-07-09 Run 2	07-Mar-200
11	Barcode:							3-9	07-Mar-200
								test template	04-Mar-200
14								test reinjections	03-Mar-200
				<b>=</b>					
-	Consumables Inf	ormation							Refres
-	Consumable	Name	Status	Days on Instrument	Expiration Date	Lot Number	Part Number		
	Polymer	POP4	384 Samples Remaining	66	01-Jan-2010 11:	51A007	4315930		
L .	Anode Buffer	AB 3xxx Buffer	299 Days Remaining	66	01-Jan-2010 02:		7275931		
	Cathode Buffer	AB 3xxx Buffer	299 Days Remaining	66	01-Jan-2010 02:		CB-431A-01		
	Capillary Array	36cm - 24 cap	105 Injections Remaining	66	01-Jan-2010 11:	80K005	4319899 - Serial # 80K2450		
•	Spatial	mation - Capillary	Array: 80K2450 Calbration Date: 03-Mar-20	09 02:53:38 PM					
	Spectral								
1	Dye Set	Chemistry Standard	Calibration	Date	Run ID				

#### Figure 6 Load Plates for Run

2. Review the consumables information and the calibration information and ensure the status is acceptable for a run.

 Enter a Run Name or use the default run name: <Start Instrument Run Date/Time Stamp> YYYY-MM-DD-hh-mm-ss-SSS (milliseconds), for example, "Run 2009-02-05-15-03-42-096" where the run start date is February 5 2009 and the run start time is 15:03:42:096.

**Note:** An instrument run begins when you click Start Run (on the Load Plates for Run screen) and ends when the last injection on the last plate has completed. For example, if you link two plates, then start the run, both plates and any duplicate injections or re-injections are part of the same instrument run. An injection is an instance of 8 or 24 samples (depending on instrument configuration) processed simultaneously under the same conditions.

**IMPORTANT!** It takes, approximately, 10 seconds for the instrument to initialize after the instrument door is closed. Do not start a run until the instrument status light is green.

When you access the Load Plates for Run screen by clicking Load Plates for Run on the Assign Plate Contents screen, the plate is automatically linked (indicated by the active Unlink button).

🧾 Plate Name:			
AB Applied Biosystems	Run Information You can edit the Run Name by entering text.		
Setup	* Run Name: Run 2009-03-04-15-33-16-629		Con
Define Plate Properties Assign Plate Contents	Plates on Instrument		
🌉 Run Instrument	Plate A (96 wells)	Link Plate	(Unlink) P
Load Plates for Run	Name: Plate 01		
Preview Run	Type: HID		
Monitor Run	Barcode:		
Review Results			

- 4. If needed, click **Unlink**, then follow the steps in "If a plate is not linked" below.
- 5. As needed, click **Switch Plates** () to assign the plate to the other position in the autosampler.
- 6. Click either of the following:
  - Create Injection List Displays the Preview Run screen where you can modify the injection list before starting the run. Go to "Review and modify the injection list in Preview Run" on page 59.
  - Start Run Displays the Monitor Run screen. Go to "Monitor the run" on page 61.

If a plate is not linked If you access the Load Plates for Run screen from the navigation pane, a plate may not be linked (indicated by the active Link button).

🛄 Plate Name:		
AB Applied Biosystems	Run Information You can edit the Run Name by entering text.	
Setup	* Run Name: Run 2009-03-04-15-33-16-629	Con
Define Plate Properties	Distance Testimores	
Assign Plate Contents	Plates on Instrument	$\frown$
🌉 Run Instrument	Plate A	Link Plate Unlink P
Load Plates for Run		
Preview Run		
Monitor Run		
Review Results		
View Conversion Results		

To link a plate:

1. Click Link Plate to display the Select Plate from Library dialog box.

L	S	elect Plate From Library								
Ir	Instructions									
2	Select row from table and click on "Link Plate" button.									
	F	ilter: Sequencing	💙 S	earch: All						
		Plate Name	T	Description						
		Place Name	Туре	Description						
	1	3-07-09 Run 1	Sequencing							
	2	3-07-09 Run 2	Sequencing		~					
	<				>					
ſ		Close		Link F	lata					
U		Close			nace					

- 2. Select a plate, then click Link Plate.
- 3. Do either of the following:
  - Click **Create Injection List**, then go to "Review and modify the injection list in Preview Run" on page 59.

or

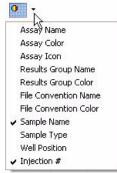
• Click Start Run, then go to "Monitor the run" on page 61.

### Review and modify the injection list in Preview Run

The Preview Run screen allows you to modify the injection list before you start the run.

- 1. Access the Preview Run screen (Figure 7 on page 59) from:
  - The Load Plates for Run screen by clicking Create Injection List.
  - The navigation pane by selecting **Preview Run**.
  - The Dashboard by clicking the **Main workflow arrow (**, then selecting **Preview Run** in the navigation pane.
- 2. Click the icon above the plate to specify the attributes to display in the plate view.
- 3. Click the plate tabs to display Plate A or Plate B.





	4 injections created -	4 in Plate A - 0 in Pla	ste B		
3	Type Assay		Instrument Protocol	Plate	Plate A Plate B 🔍 🕫 😥
0	1 IF+Norm	_POP4_xl	HID36_POP4xl_G5	3-09 Run 2	
	2 IF+Norm	_POP4_xl	HID36_POP4xl_G5	3-09 Run 2	1 2 3 4 5 6 7 8 9 10 11 12
	3 IF+Norm	_POP4_xl	HID36_POP4xl_G5	3-09 Run 2	A sample
	4 IF+Norm	_POP4_xl	HID36_POP4xl_G5	3-09 Run 2	1 1 1 2 2 2 3 3 3 4 4 4
					B sample sa
4 —					c sample
					D sample
					sample
					E 1 1 1 2 2 2 3 3 3 4 4 4
					F sample
					annale
					G 1 1 1 2 2 2 3 3 3 4 4 4
					G     1     1     1     2     2     3     3     4     4       sample samp
					<sup>6</sup> 1 1 1 2 2 2 3 3 3 4 4 4
					G     1     1     1     2     2     3     3     4     4       sample samp
					G     1     1     1     2     2     3     3     4     4       sample samp
	Legend	ion 🔮 Re-Injection	1		G         1         1         1         2         2         2         3         3         4         4         4           H         sample
			n		G         1         1         1         2         2         2         3         3         4         4         4           H         sample
	Duplicate Inject		Satus	Days on Instrument	Image: Sample
	Consumables Inf	ormation		Days on Instrument 66	I       I       I       2       2       3       3       4       4         H       sample
	Consumables	formation Name	Status		3     1     1     2     2     3     3     4     4       sample sam
	Consumables Inf	Name POP4	Status 384 Samples Remaining	66	9     1     1     2     2     3     3     4     4       sample sampl
	Consumables Inf Consumables Inf Polymer Anode Buffer	Name POP4 AB 3xxx Buffer	Status 384 Samples Remaining 299 Days Remaining	66 66	9         1         1         2         2         3         3         4         4           sample
	Consumables Inf Consumables Inf Polymer Anode Buffer Cathode Buffer	Name POP4 AB 3xxx Buffer AB 3xxx Buffer	Status 384 Samples Remaining 299 Days Remaining 299 Days Remaining	66 66 66	9         1         1         2         2         3         3         4         4         4           sample s
	Consumables Inf Consumables Inf Polymer Anode Buffer Cathode Buffer	Name POP4 AB 3xxx Buffer AB 3xxx Buffer	Status 384 Samples Remaining 299 Days Remaining 299 Days Remaining	66 66 66	9         1         1         2         2         3         3         4         4         4           sample s
	Consumables Inf Consumables Inf Polymer Anode Buffer Cathode Buffer	Name POP4 AB 3xxx Buffer AB 3xxx Buffer	Status 384 Samples Remaining 299 Days Remaining 299 Days Remaining	66 66 66	9         1         1         2         2         3         3         4         4         4           sample s

Figure 7 Preview Run screen

The Preview Run screen contains an injection list and a plate view. The injection list is linked to the plate view. Click an injection to select the associated wells in the plate view.

**IMPORTANT!** If the injection list is blank, make sure that you clicked Create Injection List on the Load Plates for Run screen.

4. To modify the injection list at any time before a run or during a run, select an injection, then click **Move Up**, **Wove Down**, and **Delete** as needed.

**Note:** Samples with assays that specify more than one instrument protocol are listed one time in the injection list for each instrument protocol.

5. To specify a duplicate injection (a replicate injection that uses the same instrument protocol as the original injection), select an injection, then click 🕎.

Sample data files for each duplicate injection can be saved in a separate folder in the results group folder if specified in the results group. For more information, see "Results group example 3: store re-injections in separate folders" on page 162.

**Note:** To use a different protocol for a replicate injection, specify a re-injection in the Monitor Run screen after you start the run.

### Start the run

When the injection list is configured, click **Start Run**. The Monitor Run screen is automatically displayed.

**IMPORTANT!** You must specify re-injections before the run completes.

**Note:** It takes, approximately, 10 seconds for the instrument to initialize after the instrument door is closed. Do not start a run until the instrument status light is green.

### Monitor the run

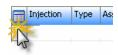
The Monitor Run screen (Figure 8 on page 61) is automatically displayed when you click Start Run in the Load Plates for Run screen or the Preview Run screen. The current injection is highlighted in green in the plate view. The injection list is linked to the plate view. Click an injection to select the associated wells in the plate view. A selected injection is highlighted in yellow in the plate view.

Conni	Run Na		nnected n 2009-03-08-13-28-46-90		User Name: Administrator Run Status: Running															
	<b>jection L</b> i		ails 4 in Plate A - 0 in Plate B																	
	Injection	Туре	Assay	Instrument Protocol	Plate Ar	Plate	A	Plate I	в							0	-		10 F	•*•
1			IF+Norm_POP4_xl	HID36_POP4xl_G5	3-09 Run 2				-			_				_				-
2			IF+Norm_POP4_xl	HID36_POP4xl_G5	3-09 Run 2	Г	1	2	:	3	4	5	6	7	8	9	10	11	12	1
3			IF+Norm_POP4_xl	HID36_POP4xl_G5	3-09 Run 2	A	4 5	ample s	ample	sample	sample	sample			samp	le sample	samp	le sampl	e sampl	le
4			IF+Norm_POP4_xl	HID36_POP4xl_G5	3-09 Run 2	ΗĤ		ample c	1 accole	1 cample	2 cample	2 cample	2	3	3	3 le sample	4	4 Is camel	4	
						E	3	ampers L 1	anipic 1	1	2	2	2	3	3	3	4	4	4	ic i
								angle o	-	-	2	2	2	3	3	3	. samp	le sampl	ie sampl	ie
								ample s	ample	sample	-	-	-	~	~	le sample	e samp	le sampl	le samp	le
						Ľ	<u>'</u>		1	1	2	2	2	3	3	3	4	4	4	
						E	5	ample s	ample 1	sample 1	2 sample	2 sample	2 sample	sample 3	samp 3	k sample 3	samp 4	le sampl 4	e sampl 4	æ
						F	5	ample s	ample	sample	sample	sample				le sample	samp	le sampl	ie samp	le
						ΗĤ		i i	1 neucla	1	2 comple	2	2	3	3	3 le sample	4	4	4	
						0	5 <mark>5</mark>	ampie s	anpe 1	1	2	2	2	3	3	3	4	4	4	c
						ŀ	1 1	ample s	ample 1	sample 1	sample 2	sample 2		sample 3	samp 3	le sample 3	samp 4	le sampl 4	e sampl 4	ie
<					>	- II -	_													



**Note:** Samples with assays that specify more than one instrument protocol are listed one time in the injection list for each instrument protocol.

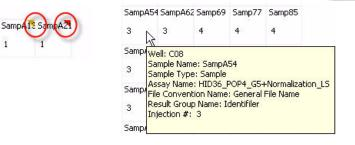
1. Click the Table Settings button, then specify the columns to show or hide in the injection list.



- 2. Optional:
  - Click the icon above the plate to specify the attributes to display in the plate view. In addition to the attributes available in Preview Run, a Flag attribute is available.

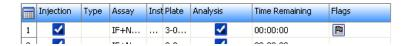
If you select the Flags attribute, yellow or red marks are displayed for wells with an Average QV value (sequencing) or an SQ value (fragment/HID) in the Fail or Suspect range. Red marks are displayed for wells with offscale data.

• Place the mouse pointer over a well to display sample details.



# Check sequence or sample quality and specify re-injections

When an injection is complete, it is flagged with  $\checkmark$  in the Injection and Analysis columns. If the software detects a problem with offscale data or low quality samples, the injection is also flagged with  $\bowtie$ .



**Note:** If the Injection, Analysis, or Flags columns are not displayed, you can click the Table Settings button, then show them in the injection list.

Injection Type As:

Assay Name

Assay Color

Assay Icon Results Group Name

Sample Type

Well Position

 Injection #

✓ Flag

Results Group Color

**File Convention Name** 

File Convention Color Sample Name

### Check sequence or sample quality

1. Expand the Flag pane at the bottom right of the screen.

Run M	atus: Cor Iame: Rur	nected 2009-03-0	08-13	3-28-46	-903	User Name: Admi Run Status: Comp			Estimat		-	ime: 07 ning: 00		09 11:2	7:35 A	M	
Injection	List Deta	ails															
njections o	reated - 8	in Plate A	- 0 in	Plate E	3												
Injection	n Type	Assay	Inst	Plate	Analysis	Time Remaining	Plate A	Plat	te B				0	•		++++	ī
		IF+N		3-0	<b>V</b>	00:00:00							-				
2 🔽		IF+N		3-0	~	00:00:00		1 2	2 3	4 5	5 6	7	8 9	10	11	12	
3 🖌		IF+N		3-0	-	00:00:00	A		samp sam								
•		IF+N		3-0	-	00:00:00			1 1 same sam		2 2 same sa		3, 5 3				
5 🗹	<u>₽₽</u>	IF+N		3-0	✓	00:00:00	В	1	1 1	2	2 2	3, 5	3,5 3	, 5 4, 6	5 4, 6	4,6	
· 🔽	2	IF+N		3-0	✓	00:00:00	C		samp sam		samp sa 2 2		same s				
Image: Second	₩	IF+N		3-0	✓	00:00:00	D	samt s	samp sam		samt sa	mp samp		amp sam	¢ sam¢	samp	
	2	IF+N		3-0	✓	00:00:00	E	samp s	samp sam	¢ sam¢ :	samp sa	mp samp	same sa	mc sam	¢ sam¢	samp	
							II F	1 samr 4	1 1 same sam		2 2 camr 42		3,5 3				
							F	1	1 1	2	2 2	3, 5	3, 5 3	, 5 4, 6	5 4, 6	4,6	
							G	samt s	samp sam		samt sa 2 2		samp si 3, 5 3				
							н		samp sam			mp samp	samp sa	amp sam	¢ sam¢	samp	
							11 -		1 1	2	2 2	3, 5	3, 5 3	, 5 4, 0	0 4,0	4,0	
Legend							Nar	a 2.00	Run 2			Barcoo	la ·				
				d 🔀 A	borted 🛃 Co	mpleted 👱	real	ie: 3*0:	PROFILE			barcoc	ю.				
te-Injection	n 🔛 Dupli	cate 📳 Fl	lags														_
Instrume	nt Run V	iews and	Flag	ç												K	Ē
inser unic	inc reality	ich's and	Thory	-											E	16	í
ray Sampl	e EPT												85	*		ď	1
												100				₽.	_
jection: 3		100	5		1200	1000 200			200		_						ğ
2			_		(seconds)				200	~ <u>Co</u>	lor	Name					ŝ
				apood (	, country						۲	EpVol	tage(kV	)			Flans Found
									>		-						n,
(																	~

The flag table displays a quick preview of sample quality and identifies samples that may need investigation.

The flag table is linked to the plate view. Click a flag to select the associated well in the plate view:

**Note:** If no samples are listed in this pane, no flags were found and the samples have passed quality checks.

• All samples passed

sample sample sample sample 3, 5, 7, 13, 5, 7 3, 5, 7, 4,6 2 Flags Found Display All Flags ~ 0 InjecOffscale SQ/QV Cap # Samp 3 **F** 9 samp

v 🕜

9

18

sample

sample

.....

InjecOffscale SQ/QV Cap # Sample N

2 Flags Found

Display All Flags

3

3

0, 0

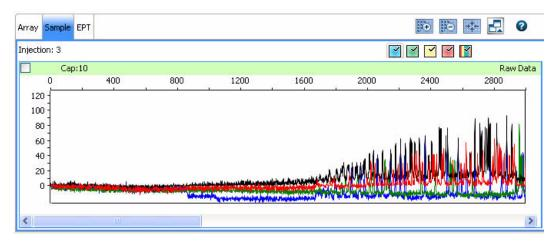
- E At least one sample is in the suspect range and requires review
- At least one sample is offscale or is in the suspect range

2. To filter the flag table, select a flag type. To display HID flags, select **All**. To sort the table, double-click column headers.

The flags you may see in the flag table are:

Flag/Symbols	Description
Offscale	[2] (red) At least one data point in the analysis range has saturated the CCD camera.
(green or red)	Note: In the View Results screen, an offscale sample is flagged with $\underline{\mathbb{A}}$ .
Average Quality Value (sequencing)	(yellow) or [] (red) The Average Quality Value (based on CRL, Trace Score, and QV20+ results) is in the Suspect or Fail range. For information, see "Basecalling protocol – QV settings" on page 178.
(green, yellow, red)	
Sizing Quality (fragment/HID)	(yellow) or [1] (red) The Sizing Quality is in the Suspect or Fail range. For information, see, Table 15 on page 183 or Table 17 on page 188.
(green, yellow, red)	<b>IMPORTANT!</b> Normalization is not applied to samples with (red) Sizing Quality.

3. Click a row in the flag table, then click the Sample tab in Instrument Run Views to display the associated data in the Sample view.



#### Specify re-injections

You can specify a re-injection before the run completes. A re-injection physically re-injects all samples in the capillary array. You can select a different instrument protocol than the original injection and can specify whether to collect data for all or only selected samples in the array.

1. Select the injections or wells to re-inject:

**Note:** Re-inject is grayed if you select an injection that contains more than one results group, or if you select flags in the flags table that correspond to samples with different results groups. To enable Re-inject, select samples that specify the same results group.

To collect data for all wells in an injection	1. Select the injection in the injection list.
	2. Click 🕎 Re-inject.
To collect data for only specific wells	1. Select the injection.
(Samples with assays that specify more than one instrument protocol are listed one time in the injection list for each instrument protocol)	2. Select in the array view the capillary that corresponds to the well or sample of interest (see "Array view" on page 77).
<b>Note:</b> You can also specify re-injections for specific samples in Review Results.	3. Click 🕎 Re-inject.
To collect data for only samples that contain flags	1. Select the samples in the flag table (see "Check sequence or sample quality" on page 62).
	2. Click 🕎 Re-inject.

**Note:** If you are running an HID plate, see "Re-injections of HID allelic ladder samples" on page 67.

- 2. In the Re-injection dialog box, select options, then click OK:
  - The protocol to use for the re-injection: original, modified, new, or one from the library
  - When to make the re-injection

**Note:** Sample data files for each re-injection can be saved in a separate folder in the results group folder if specified in the results group. For more information, see "Results group example 3: store re-injections in separate folders" on page 162.

U Re-injection	
Re-injection	16
Create a Re-injection	11
	Ø
Instrument Protocol Options	
• Reuse the existing protocol	
O Modify the existing protocol	
O Create a new protocol using the template:	HID36_POP4xl 🔽
OReuse a protocol in the library:	HID36_POP4xI_F
Placement of Re-injections	
<ul> <li>Following all injections</li> </ul>	
O After original injection	
	OK Cancel

#### If you select a protocol other than the original

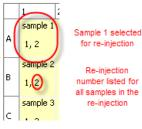
If you select a protocol other than the original, the software:

- Creates a copy of the assay specified for the re-injected well (Original\_Assay-1).
- Adds the new or modified instrument protocol to Original\_Assay-1.
- Assigns Original\_Assay-1 to the re-injected well only. ٠
- Saves the plate (the software does not save the copy of the assay to the library). ٠

#### How re-injections are displayed in the plate view

If the Injection Number attribute is selected for display in the plate view, the number of the original injection and the re-injection are shown.

Note: If you select only specific wells for the re-injection (which physically re-injects all samples for the capillary array but collects data only for the selected wells), the re-injection number is displayed for all samples in the re-injection, not just the samples selected for data collection.



#### Re-injections of HID allelic ladder samples

If you select to re-inject a sample that includes an allelic ladder in its results group, but the allelic ladder is not part of the injection, the software prompts you to select one or more allelic ladder samples to re-inject.

For example:

- You are running an 8-capillary instrument, and you have specified one results group for each set of three injections (for more information, see "Results group example 2: store one allelic ladder per run folder (8-capillary instruments)" on page 161)
- The allelic ladder sample is in Injection 1.
- You select for re-injection a sample that is in injection 2.
- The software prompts you to select one or more allelic ladder samples to re-inject.

The allelic ladders available to select are from the same plate and within the same results group as the original injection. If the results group does not contain an allelic ladder sample, the software does not prompt you to select one for re-injection.

Je Add Allelic Ladder to Re-injection	×
Add Allelic Ladder to Re-injection	-
	0
Select zero (0) or more Allelic Ladder samples.	
A01_allelic ladder	
Deselect All	
<ul> <li>Allelic Ladder Options</li> <li>Collect data for all other wells in the selected allelic ladder's injection</li> <li>Apply modified instrument protocol to all wells in the selected allelic ladder's injectio</li> </ul>	n
OK Cancel	

In the Add Allelic Ladder to Re-injection dialog box:

1. Select one or more allelic ladder samples.

**IMPORTANT!** The software does not display the well location of allelic ladder samples in this dialog box. To identify allelic ladder samples for re-injection, include the well position in the allelic ladder sample name when you assign plate contents.

2. Select whether to collect data for the remaining samples in the allelic ladder re-injection.

3. Select whether to apply a modified instrument protocol to the allelic ladder re-injections, or whether to use the original instrument protocol for the allelic ladder re-injection(s). You will select the modified protocol in the next screen.

**IMPORTANT!** Allelic ladders that are injected under the same conditions are recommended to accurately genotype samples in the secondary analysis software (GeneMapper<sup>®</sup> *ID-X* Software v1.2 or later).

- 4. Click OK.
- 5. Specify the remaining re-injection settings as described in "Specify re-injections" on page 65.

Two re-injections are added to the injection list. The first re-injection collects data for the selected sample. The second re-injection collects data for the allelic ladder.

### **Review completed injections in Review Results**

You can review results for any completed injections. Select the injection, then click **Review Results**. The samples for the injection are loaded in the Samples Table in Review Results. For more information, see "Review Results" on page 79.

### Start and stop a run

Start a run	You can start a run in the:
	• Load Plates for Run screen (see "Load plates for run and create the injection list" on page 56).
	• Preview Run screen (see "Start the run" on page 61).
Pause and	As needed, click:
resume a run	• <b>D</b> Pause – Pauses the run after the current injection completes (the <b>D</b> symbol is not displayed in the injection list because the injection continues to completion).
	• <b>Resume</b> – Resumes the run.
Abort or	As needed, click:
terminate	• <b>Solution</b> Abort – Stops the current injection. Do not click Delete to stop an injection.
	<b>IMPORTANT!</b> You can stop the current injection only when the front panel indicator is blinking green. If you click Abort when the front panel indicator is solid green, the physical injection is already completed (although the software is still processing the information) and a message is displayed indicating that there is no injection in process.
	• <b>Solution</b> Terminate – Stops the instrument run. Terminate is active only when a run is paused.

## More features in Assign Plate Contents

### Use the Plate View

Name samples in	To name samples in th	e Plate View:						
the Plate View	To name one sample	<ul> <li>Click a well, then type a sample name directly into the field, then press Enter.</li> <li>or</li> <li>Copy and paste a name from another well.</li> </ul>						
		To set the directio cursor when you p • Click Row to	oress Enter: o set the En		1			
		<ul> <li>key to move th vertically to the</li> <li>Click Column Enter key to me horizontally to column.</li> </ul>	2					
	To name multiple samples	<ol> <li>Click a named</li> <li>Click-drag multication</li> <li>Right-click and selected fields</li> </ol>	tiple wells.	or <b>Fill Se</b> i	ries to popula	ate the		
		<b>Note:</b> To use Fill Series, type a number as the last character of the named well). You can also copy and paste sample names.	A Sample 1 B	2	3	4		
			C D E	Cut Copy Paste Fill		Ctrl+> Ctrl+( Ctrl+\ Ctrl+I		
			F		Fill Series Fill Sample N Delete Rename Sam			
	To name all wells at one time	<ol> <li>Select all wells</li> <li>Select assays, conventions, a group for the p</li> <li>Enter name and sample type (in Customize San pane) for the w</li> </ol>	file name nd results late. d select i the nple Info			7     8     9     10     11       1     1     1     1     1       1     1     1     1     1       1     1     1     1     1       1     1     1     1     1       1     1     1     1     1       1     1     1     1     1       1     1     1     1     1       1     1     1     1     1       1     1     1     1     1       1     1     1     1     1		

Sample7 Sample8

Customize the plate view	<ul> <li>Click Show In Wells to specify the attributes to display in wells.</li> <li>Click Select Wells to select wells with a specific attribute.</li> </ul>	<ul> <li>Show In Wells</li> <li>Assay Name</li> <li>Assay Color</li> <li>Assay Icon</li> <li>Results Group Name</li> <li>Results Group Color</li> <li>File Convention Nar</li> <li>File Convention Color</li> </ul>			
E Zoom II	Click Zoom In, Zoom     Out, and Fit as     needed.	✓ Sample Name Sample Type Well Position	Select Wells     Assay Name     Sample Type     Sample Name     Results Group Name     File Convention Name     Well Position	Array Selection Sample1 Sample2 Sample3 Sample4 Sample5 Sample6	₹73 974

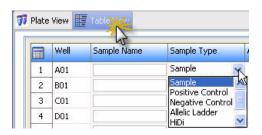
View the capillary/plate map

Click **Array Selection** to select wells by injection. Click again to turn off array selection.

t Wells 🔻	Array Sele	tion Ro	w 🔁 Colum
7	8	9	10
_			

### Use the Table View

- 1. Click Table View.
- 2. Click the Sample Name field, then type a name.
- Click next to each field, then select a setting.



4. Right-click a column header, then select **Fill** or **Fill Series** to populate the selected fields (to use Fill Series, type a number as the last character of the named well).

	Well	Sample Name	Sample Typ	be	Assa
1	A01	Sample 1	Cut	Ctrl+X	
2	B01		Copy Paste	Ctrl+C Ctrl+V	
3	C01		Fill	Ctrl+D	
4	D01		Fill Series		
5	E01		No Data	Y	_
6	F01	1	No Data		

Note: You can

double-click column

headers to sort columns. Multi-column sorting is supported (see "Multi-column sorting" below).

### Sort and customize tables

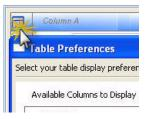
Multi-column sorting

- You can sort any table in the software. Multi-column sorting is supported:
  - Double-click a column header to sort the column.
  - Alt+Shift-click another column header to sort another column.
  - Alt+Shift-click a third column header to sort a third column.

Numbers in the column headers reflect sort order.

		~		-		0
	Column A	(-1)	Column B	( 2 )	Column C	- 3
1		$\sim$		$\mathbf{\nabla}$		$\sim$
2						

**Customize tables** You can customize any table in the software. Click the Table Settings button, then specify the columns to show or hide.



### Add assays, file name conventions, and results groups to a plate

1. If no assay is listed at the bottom of the Assign Plate Contents screen, add at least one assay. You can specify different assays for different wells.

	Assays			File	Name E
		Actions	•		
	Add from Library	<			Add fro
	Create New Assay	2		Create I	New File
A	dd Assay From Library				
	ct rows from table and click on				
	ct rows from table and click on ilter: Fragment	"Add To Plate"	button. Search:	All	
			Search:	All Internet Protocol	Prima
	ilter: Fragment		Search:		
F	ilter: Fragment Assay Name	▼ Type Fragment	Search: Instru Fragr	Iment Protocol	1 Fragn

2. (Optional) If no file name conventions or results groups are listed at the bottom of the Assign Plate Contents screen, add as needed. File name conventions and results groups are optional, but are very useful for naming and organizing data files.

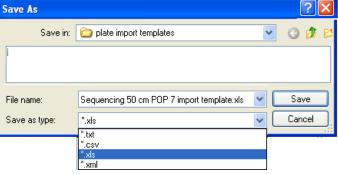
### Create a plate for importing

Create a plate import template	The 3500 Series Data Collection Software allows you to import plate information from files that you create in an application other than the 3500 Series Data Collection Software.
	To create a template for importing plate information, set up a plate in the 3500 Series Data Collection Software, then export it to create a file that contains the correct header and column information for importing:
	1. In the Dashboard, click <b>Create Plate from Template</b> .
	2. In the Open Plate Template from Library dialog box:
	a. Select a filter to display the plate template type of interest.
	b. Select a plate template, then click <b>Open</b> .
	3. Enter a name for the plate, then specify the capillary length and polymer type for the plate.
	4. Click Assign Plate Contents.

5. In the Assign Plate Contents screen, click 🛃 Export.

**Note:** Before you click Export, you can assign other plate elements to the plate import template as described in "Assign plate contents" on page 46.

- 6. Select a file type for the plate import template.
- 7. Enter a name and location for the plate record template.



8. Click Save.

The figure below shows the format of the exported plate.

	A	В	С	D	E	F	G	Н	
1	3500 Plate Layout Fil	e Version 1.0							
2									
3	Plate Name	Application Type	Capillary Length (cm)	Polymer	Number of Wells	Owner Name	Barcode Number	Comments	
	plate import template	Sequencing	50	POP7	96				
5									
6	Well	Sample Name	Assay	Results Group	File Name Convention	Sample Type	User Defined Field 1	User Defined Field 2	User Defi
7	A01								
8	B01								
9	C01								
10	D01								

# Create a plate import file

- 1. Open a plate import template (see "Create a plate import template" on page 73).
- 2. Save the plate import template under a new name.
- 3. Enter sample names (required).
- 4. (Optional) Enter information for the remaining columns.

**Note:** If you specify assay, results group, or file name convention names, the names you enter must exactly match the names of existing items in the library.

5. Save the plate import file.

#### Edit a plate

You can edit a plate from:

- Library Select a plate, then click **Z** Edit.
- Dashboard Click Edit Existing Plate.
- Define Plate Properties screen Select Open Plate > Edit Existing Plate.
- Assign Plate Contents screen Select Open Plate > Edit Existing Plate.

### Import and export a plate

You can import and export plates from:

- **Plates library** Plates in .xml format for use on another 3500 or 3500xL analyzer instrument. See "Import and export a library entry" on page 141.
- **Define plate properties** Plates in .txt, .csv, and .xls format files you create that contain plate information in a specific format.
- Assign Plate Contents Plates in .txt, .csv, and .xls format files you create that contain plate information in a specific format.

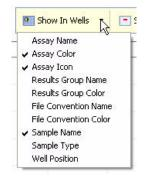
#### Create a plate template

A plate template contains default settings that you can edit when you create a plate from the template.

- 1. Create a plate (see "Create a new plate" on page 144).
- 2. (Optional) Add sample names and sample types (see "Name samples and assign sample types in the plate view" on page 48).
- 3. (Optional) Add the assays, file name conventions, and results groups appropriate for this plate template's application (see "Add assays, file name conventions, and results groups to a plate" on page 73).

Adding assays, file name conventions, and results groups to the plate template automatically displays these items in the Assign Plate Contents screen when you open the plate template. You do not have to add these items from the library for each plate you create.

- (Optional) Click **Show In Wells** to specify the attributes to display in wells in the template.
- Select Save Plate ➤ Save As Template. The software displays the template icon below the plate layout.



### Specify the default plate type for the Open Plate dialog box

Library Maintenance Tools \* Manage \* Preferences Help \* Log Out 2 🔚 Preferences type filter text not used **Plate Setup** 🖃 System Choose the default application type, polymer type, and capillary length to use when the ins Date Format Instrument Settings Application Sequencing ✓ Polymer Type POP7 ✓ Capillary Length 50 Scheduler Preference 😑 Sequencing Settings Export Spectral Calibration 🖃 User Plate Setup **Reports Settings** Run Setup Sequencing Settings Trace Trace Print Choose the default Assign Plate Contents View. Trace Quality Trace Quality Reports Assign Plate Contents View Table Plate h Table

Specify the default plate type for the Open Plate dialog box in Preferences.

### Save electronic version of reports

When you print any report, you can select **CutePDF Writer** as the printer, to save the report to .pdf.

### More features in Load Plate for Run

#### Link a plate from the recent plates or recent runs tab

Instead of clicking Link to select a plate, you can click-drag a plate from the Recent Plates tab (pending plates) or the Recent Runs tab (processed plates).

Plate Unlink	Plate B	Link Plate	Unlink	Recent Plates	Recent Runs
				Name	Date Modified
				3-07-09 Run 2	0

### More features in Monitor Run

### **Review the Instrument Run views**

Select an injection, then click an instrument run view tab. As needed:

- Click 😥 🔁 🐝 to zoom in and out
- Click I to detach a view and display it in a separate window that you can move around on the screen.

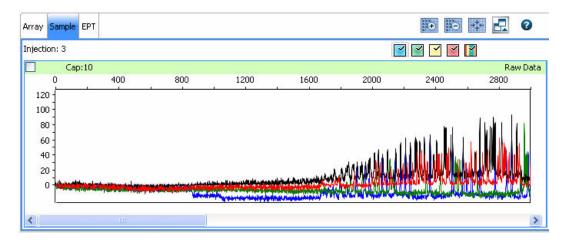


To locate a detached view, click the 3500 task bar icon.

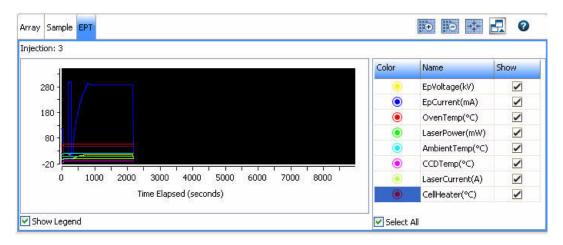
**Array view** The Array view shows the color data (based on the dominant fluorescence color) for each capillary as a function of instrument scan number (time). Adjust the brightness and color by using the slider bars above the view.



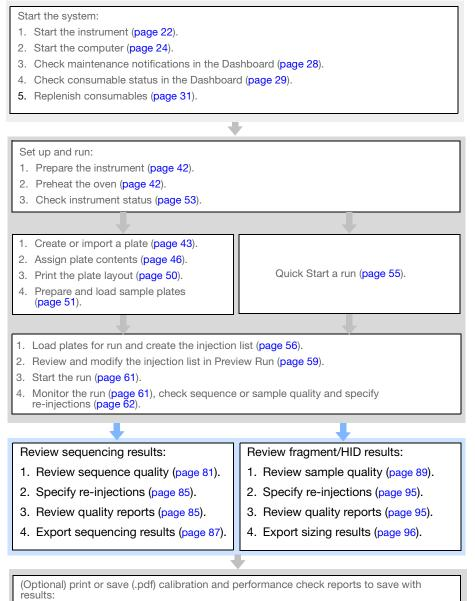
**Sample view** The Sample view shows the relative dye concentrations as a function of instrument scan number (time) for the selected capillary. You can select and deselect the dye colors to display.



**EPT view** The EPT view (ElectroPhoresis Telemetry) shows instrument data conditions (laser power, temperatures, electrophoresis voltage) as a function of time. In the legend to the right of the EPT view, you can select and deselect the traces to display in the view.



### Workflow



- Spatial calibration (page 99)
- Spectral calibration (page 103).
- Sequencing install standard performance check (page 122).
- Fragment or HID install standard performance check (page 132).

### **Review Sequencing Results**

### Access the View Sequencing Results screen

Access the View Sequencing Results screen from:

- The Monitor Run screen by clicking **Review Results**.
- The navigation pane by selecting View Sequencing Results.
- The Dashboard by clicking View Run Results.





#### Review results for the currently running plate

If you access the View Sequencing Results screen while an instrument run is in progress, the Trace Quality View lists results for completed injections in the current run.

Select one or more samples, then click by Open Trace to display their data in the Trace pane.

**Note:** The basecaller version listed in the basecalling protocol is limited to a 3-digit number. The version listed in sequencing results is a 4-digit number. The fourth digit is an internal number used by the software.

	ace Quality View	≚ Trac	e Quality Rep	orts Sele	ct Table:	Analysis Result	5 💌	Select F	Filter: A	II Traces	~	
												i
	Sample Filename		Assay Name		PA Prot	ocol Name	Basecall Version	n 🛛	Mobility	File	В	ase Sp 🔨
1	B04_hCV111684	29	StdSeq50_PC	DP6_v3.1	BDTv3.	1 PA Protocol	KB 1.4.1.6		KB_3500	_POP6_BD	T 1	3.81
2	B09_hCV111923	13_9	StdSeq50_PC	DP6_v3.1	BDTv3.	1 PA Protocol	KB 1.4.1.6		KB_3500	_POP6_BD	T 1	3.76
3	C01BAC7_v3.1ft	wd7	A2_v3.1Rapi	idSeq50	v3.1_S	equencing	KB 1.4.1.4		KB_3500	_POP7_BD	T 1	4.18
4	C05BAC2_v3.1ft	wd8	A1_v3.1Shor	tReadS	v3.1_5	equencing	KB 1.4.1.4		KB_3500	_POP7_BD	T 1	4.01
5	B04_hCV111684	29	StdSeq50_PC	DP6_v3.1	BDTv3.	1 PA Protocol	KB 1.4.1.6		KB_3500	_POP6_BD	T 1	3.81
6	B09_hCV111923	13_9	StdSeq50_PC	DP6_v3.1	BDTv3.	1 PA Protocol	KB 1.4.1.6		KB_3500	_POP6_BD	T 1	3.76
7	C01BAC7_v3.1ft	wd7	A2_v3.1Rapi	idSeq50	v3.1_S	equencing	KB 1.4.1.4		KB_3500	_POP7_BD	T 1	4.18
8	C05BAC2_v3.1ft	wd8	A1_v3.1Shor	tReadS	v3.1_5	equencing	KB 1.4.1.4		KB_3500	_POP7_BD	T 1	4.01 🖉
<	1											>
												_
🖻 Tra	асе		- 💹	AA   <u>E</u>		ACG	1 1 1	Sel	: Tab Ke	y to:	-   3	) 🗩 j
6 BO4	4 hCV11168429 :	1 🔀	B Copy (1	.0) of B09		ab Copy (10) of	C01B a	Copy ()	10) of C	D58	» <sub>8</sub>	
	4_hCV11168429_		💩 Copy (1	.0) of B09		💩 Copy (10) ol	C01B a	🖻 Copy (	10) of C	D5B	»,	
	inates (x,y): 1	151,9	]								•	
		151,9	Copy (1			ab Copy (10) of 50 2000 205		Copy (:	10) of C 2250	2300 23	•	-
	inates (x,y): 1 1650	151,9 1700	  750 1800	1850	1900 19	50 2000 205	0 2100 2150	2200			•	- (
	inates (x,y): 1 1650	151,9 1700	]	1850	1900 19	50 2000 205	0 2100 2150	2200	2250	2300 23	•	- (
	inates (x,y): 1 1650	151,9 1700	] 1750 1800 T CCAA GG0	1850	1900 19	50 2000 205	0 2100 2150	2200 	2250	2300 23	50	- (
Coord	inates (x,y): 1 1650 50 CTTTGTG	151,9 1700 GACT G	] 1750 1800 T CCAA GG0	1850 JULIUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUU	1900 19	50 2000 200 NTTT GT GGC/	0 2100 2150	2200 	2250 	2300 23	50	- (
Coord	inates (x,y): 1 1650 50 CTTTGTG 140	151,9 1700 GACT G	] 1750 1800 T CCAA GG0	1850 JULIUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUU	1900 19	50 2000 200 NTTT GT GGC/	0 2100 2150	2200 	2250 	2300 23	50	- (
Coord	inates (x,y): 1 1650 50 CTTTGTG 140	151,9 1700 GACT G	] 1750 1800 T CCAA GG0	1850 JULIUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUU	1900 19	50 2000 200 NTTT GT GGC/	0 2100 2150	2200 	2250 	2300 23	50	- (
Coord 15	inates (x,y): 1 1650 50 CTTTGTG 140	151,9 1700 GACT G	] 1750 1800 T CCAA GG0	1850 JULIUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUU	1900 19	50 2000 200 NTTT GT GGC/	0 2100 2150	2200 	2250 	2300 23	50	- (
Coord 15	inates (x,y):	151,9 1700 GACT G	] 1750 1800 T CCAA GG0	1850 JULIUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUU	1900 19	50 2000 200 NTTT GT GGC/	0 2100 2150	2200 	2250 	2300 23	50	- (
Coord 15	inates (x,y):	151,9 1700 GACT G	] 1750 1800 T CCAA GG0	1850 JULIUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUU	1900 19	50 2000 200 NTTT GT GGC/	0 2100 2150	2200 	2250 	2300 23	50	- (
Coord 15 10	inates (x,y):	151,9 1700 GACT G	] 1750 1800 T CCAA GG0	1850 JULIUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUU	1900 19	50 2000 200 NTTT GT GGC/	0 2100 2150	2200 	2250 	2300 23	50	- (
Coord 15 10	inates (x,y): 1 1650 50 1000 CTTT GT G 140 100 1000 1	151,9 1700 GACT G	] 1750 1800 T CCAA GG0	1850 JULIUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUU	1900 19	50 2000 200 NTTT GT GGC/	0 2100 2150	2200 	2250 	2300 23	50	- (
Coord 15 10	inates (x,y): 1 1650 50 1000 CTTT GT G 140 100 1000 1	151,9 1700 GACT G	] 1750 1800 T CCAA GG0	1850 JULIUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUU	1900 19	50 2000 200 NTTT GT GGC/	0 2100 2150	2200 	2250 	2300 23	50	- (
Coord 15 10	inates (x,y): 1 1650 50 1000 CTTT GT G 140 100 1000 1	151,9 1700 GACT G	] 1750 1800 T CCAA GG0	1850 JULIUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUUU	1900 19	50 2000 200 NTTT GT GGC/	0 2100 2150	2200 	2250 	2300 23	50	- (

#### Review previously run samples

If you access the View Sequencing Results screen when no run is in progress and no plate is linked, no samples are listed. (If the plate from the most recent run is linked, the results from that plate are displayed.)

To view results for samples other than those from the most recent run, click **import**, then select the samples to review.

### **Review sequence quality**

1. Display Metric Analysis results to review sample basecalling and trimming results.

ab Trace Quality View		
Select Table:	Metrics Analysis Results	Select Filte
	Analysis Results Run Information Data Collection Information	M
	Metrics Analysis Results Trace Identification	A atocol Nar

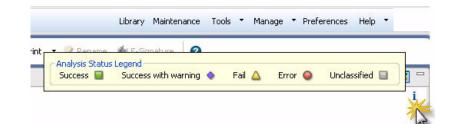
- 2. Click the Table Settings button, then specify the columns to show or hide.
- 3. Double-click column headers to sort columns. Multi-column sorting is supported (see "Sort" on page 97).



4. Review the results:

Result	Description
Trace Score	The average basecall quality value (QV) of bases in the clear range sequence of a trace.
	The clear range is the region of the sequence that remains after excluding the low-quality or error-prone sequence at the 5' and 3' ends. The clear range is calculated by the KB basecaller using QVs.
CRL	The longest uninterrupted segment of bases with a Quality Value (QV) $\ge$ 20. In addition to evaluating the QV of a base call, the software considers the QV of adjacent bases within +/- 20 bases, before including a base in the continuous read length.
QV20+	The total number of bases in the entire trace that have basecaller quality values equal to or greater than 20.
Trace Score Quality CRL Quality QV20 Quality	Pass/fail/check determined by the settings in the Basecalling protocol QV Settings tab.
PUP Score	A measure of noise as calculated as the ratio of the fluorescence signal of the highest secondary peak to the fluorescent signal of the main called base.

- 5. Review warnings:
  - a. Scroll to the right of the Metric Analysis table to display the Warning column.
  - b. Display the Analysis Status legend.



c. Review warnings:

Result	Description			
Success	Basecalling and trimming successful.			
Success with warning	Basecalling successful, trimming not successful. Warning messages are listed in the Warning/Error Message column (default position is the last column in the table).			
🛆 Fail	Basecalling and trimming failed, no results generated.			
Error	Basecalling and trimming failed due to internal software error, no results generated.			
Unclassified	No analysis performed.			

6. (Optional) Click **Minimize** and **Restore** to collapse and expand the samples table.

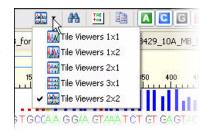
07	0
No.	Restore
Minimize	2010/01/01/01/01/01/01

### **Review traces**

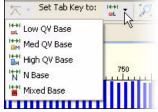
- 1. Select the samples of interest in the samples table, then click 🛅 Open Trace.
- 2. Select items from the trace toolbar to manipulate the trace as needed. Place the mouse pointer over a button for the description of the button.

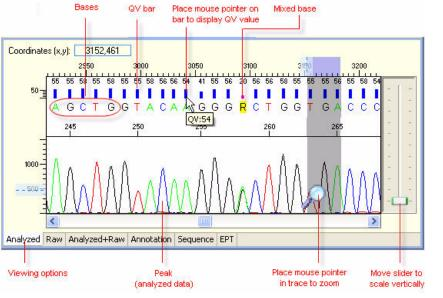


- 3. (Optional) Modify trace display:
  - Use the Tile Viewer options to display up to four traces at a time.
  - Set trace colors in Preferences (see "Set sequencing preferences" on page 36).



- 4. Set the category of base for the Tab key.
- 5. Review traces: press **Tab** to review bases from left to right in a trace. **Shift+Tab** to move right to left.





6. Click the tabs at the bottom of the trace pane for different views of the data.

2-2-2-2-2-2		~			- 2-
	<	1111	Ĵ	18	
Analyzed	Raw	Analyzed+Raw	Annotation	Sequence	EPT

### Understand Quality Values (QVs)

Quality value ranges		•		commends the ferences" on j	following ranges for QVs (set in Preferences, page 36):
	• P	ure bases	– Low	QV ≤ 15, Me	dium $QV = 15$ to 19, High $QV = 20+$ (default)
				- ·	edium $QV = 5$ to 10, High $QV > 10$ (investigate our application)
	Note:	The predic	cted pro	obability of er	From for a basecall is high $QV > 10$ .
	baseca		col (se	e "Basecallin	n (set the clear range) using quality values in the g protocols library (primary analysis –
Pure base versus mixed base QVs				ases have the ne following:	same probability of error for the associated
	• H	igh-quality	y pure	bases typically	y have QVs of 20 or higher.
		he distribu f pure base		quality value	s for mixed bases differs dramatically from that
	• Fe	or mixed b	ases, q	uality values	greater than 30 are rare.
	pi		of error	• •	ned quality values as low as 5, because the bases is higher. Review mixed bases with QVs
Quality values					
(QV) and probability of	QV	Ре	QV	Ре	
error (Pe)	1	79.0%	30	0.10%	
	5	32.0%	35	0.032%	
	10	10.0%	40	0.010%	
	15	3.2%	45	0.0032%	
	20	1.0%	50	0.0010%	
	25	0.32%	60	0.00010%	

**Display** 1. Click View Thumbnails to display results as thumbnails. **thumbnails** 

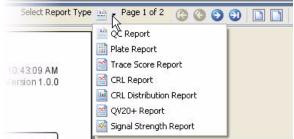
elect Filter: All Traces			
		Melhecka Willer	and a straight and and a straight and a straight and a straight a straight a straight a straight a straight a st
A04_A04.ab1	A05_A05.ab1	A06_A06.ab1	B04_B04.ab1
	all little and all allowed	Internet Hills	West and the standard and a
C06_C06.ab1	D04_D04.ab1	D05_D05.ab1	D06_D06.ab1
While and Weller while	and any filled and	talandy Mildle	
F05_F05.ab1	F06_F06.ab1	G04_G04.ab1	G05_G05.ab1
2. Sort as neede	d.	ab	Trace Quality View
-	ignal across all sa J <b>niform Y Scalin</b>	-	ort By Trace File Name
4. Click View Tapane.	ables to close the	thumbnail	Average Raw Signal Intensity Trace Score CRL Mobility File
Select Filter: A	Il Traces 🛛 💟	s:	Run Module Name 2. Run Name 2. Instrument Name Well ID

### Specify re-injections

Before the run is complete, you can select a sample, then click **M** Re-inject.

### View, print, and save (.pdf) trace quality reports

View Trace Reports 1. Click View Trace Reports to see the available reports for traces and print the reports you want. You can set defaults for the reports in Preferences (see "Set sequencing preferences" on page 36).



Well ID Capillary Results Group

2. Select the report type and review the content of each report. See "Report options" on page 86.

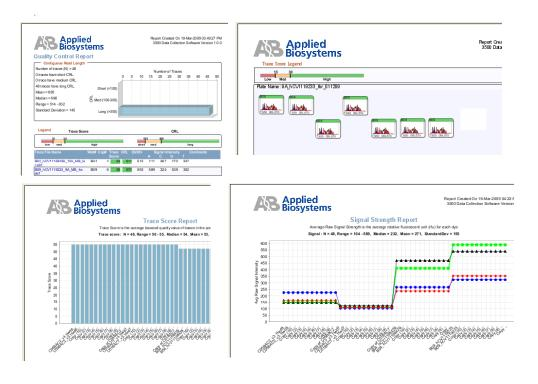
3. Modify report settings as needed. You can specify additional report settings in Preferences (see "Trace Print (user preference)" on page 38, "Trace Quality (user preference)" on page 38, and "Trace Quality Report (user preference)" on page 39).

Select Report Type 🔛 🗸 Page 1 of 2 🛛 🕝 🧿 🕥 🗋 📑 🔖 🗙 🖓
Modify report settings
Sort data based on
This setting applies to the Trace Score Report, CRL Report, QV20+ Report, and Signal Strength Report.
Run Name
Capillary Number
Signal based on
This setting applies to the QC Report, and Signal Strength Report.
Average Raw Signal Intensity
O Average Raw Signal to Noise Ratio
Font settings
Select the font to be used in reports.
Arial
10 🐱
OK Cancel

- 4. Double-click different elements in the report to open the Trace view and display the associated sample.
- 5. To print the report, click 🍓 **Print**, then preview or print.
- 6. To save the report electronically (.pdf), print the report and select **CutePDF Writer** as the printer.

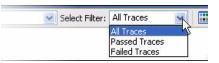


- QC One-page bar chart that shows trace score statistics and results for each selected sample.
  - Plate One-page per plate for all selected samples that shows the well-location thumbnail raw data traces with color-coded headers that reflect Trace Score quality.
  - Trace Score, CRL, and QV20+ One-page bar chart that shows trace score, CRL, or QV20+ statistics and results for each selected sample.
  - **CRL Distribution** One-page bar chart that shows CRL statistics and CRL results distribution for all selected samples.
  - **Signal Strength** One-page graph that shows with average sequencing dye signal strength for all selected samples.



### Export sequencing results

- 1. Filter the table of interest.
- 2. Select an export option: Results, Reports, or Traces.



3. Select the export options and the location for the export file, then click **OK**.

The file(s) are exported to the specified location with the following naming conventions:

- **Results** *export\_ReportName*.txt
- **Reports** *ReportName*.\* (\* is the format you selected: .txt, .xls, .pdf, .html)
- **Traces** *FileName*.\* (\* is the export format you selected: .annotation.txt, .phd.1, .scf, .fsta, .qual, .seq)

### **Review Fragment/HID Analysis results**

### Access the View Fragment/HID Results screen

Access the View Fragment/HID Results screen from:

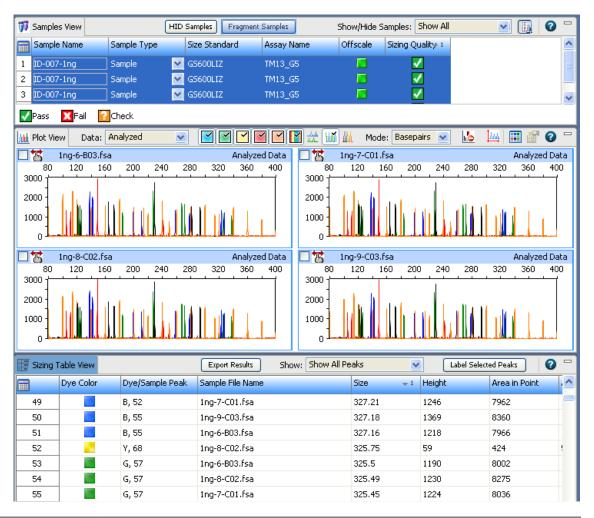
- The Monitor Run screen by clicking **Review Results**.
- The navigation pane by selecting View Sequencing Results.
- The Dashboard by clicking View Run Results.



#### Review results for the currently running plate

If you access the View Fragment/HID Results screen while an instrument run is in progress, the samples table lists results for completed injections in the current run.

Select one or more samples in the samples table to display their data in the plot view and sizing table view.



#### Review previously run samples

If you access the View Fragment/HID Results screen when no run is in progress and no plate is linked, no samples are listed. (If the plate from the most recent run is linked, the results from that plate are displayed.)

To view results for samples other than those from the most recent run, click **import**, then select the samples to review.

<b>Note:</b> By default, the Fragment Samples view is selected. If you	arts 🝷 🌺	Print 🔻 i	Import	🔊 Remove	MM Re-Inject	- Save
are importing HID files, click		HID Sam	ples	Fragment Sampl	es	Show/Hide
HID Samples.	nple Type	Size St.	ard	Assay Name	PA Pro	otocol

### **Review sample quality**

- 1. In the samples view, click the Table Settings button, then specify the columns to show or hide.
- 2. Double-click Offscale, Pull-Up (fragment), Broad Peak (HID), and SQ columns to sort suspect and failing flags to the top of the table.



Multi-column sorting is supported (see "Sort" on page 97).



Flag/Symbols	Description				
Offscale	At least one data point in the analysis range has saturated the CCD camera.				
-	Note: In the Monitor Run screen, an offscale sample is flagged with 📃.				
Spectral Pull-Up (fragment analysis only)	At least one peak contains a pull-up peak.				
	A pull-up peak is identified when the peak height of the minor peak is $\leq X\%$ of and within $\pm Y$ data point of the major peak, where X and Y are values you specify. See Chapter 6, Manage Library Resources.				
Broad Peak (HID analysis only)	At least one peak exceeds the Broad Peak threshold.				
	Broad peaks affect Sizing Quality. See Chapter 6, Manage Library Resources.				
	<b>Note:</b> The value displayed when you place the mouse pointer over a Broad Peak flag is an internal value and does not reflect the peak width.				
Normalization Limit	Sample was collected with a normalization size standard, sample Normalization Factor is within range.				
	<ul> <li>A – Sample was collected with a normalization size standard, sample Normalization Factor is not within range.</li> </ul>				
	<ul> <li>No Data – Normalization is enabled, but Sizing Quality is </li> </ul>				
	• NO – Sample was not collected with a normalization size standard.				
	• N/A – Sample was not collected on a 3500 or 3500xL analyzer instrument.				
	For more information, see "Review normalized data" on page 90.				
	<b>Note:</b> If the Sizing Quality is <b>I</b> , normalization is not applied, even if the Normalization Factor is within the normalization range.				

Flag/Symbols

Sizing Quality



**Note:** If the Sizing Quality is applied, even if the Normalization Factor is within the normalization range.

#### Description

The Sizing Quality is in the Fail or Suspect range. Place the mouse pointer over a flag to display the Sizing Quality value for the sample. See Chapter 6, Manage Library Resources.

- 3. Click a flag in the samples table, or select samples in the samples table to display the associated data in the Plot View and Sizing Table View.
- 4. (Optional) Modify the sample view:
  - Right-click the Size Standard field to view the size standard for a sample.
  - Click **Minimize** and **Restore** to collapse and expand the samples table.



### Review normalized data

Normalization corrects for instrument, capillary, and injection variability. When specified in the primary analysis protocol, the software calculates a normalization factor for each sample. The normalization factor is used as a multiplier to adjust the peak height of the sample peaks relative to the GS600 LIZ<sup>®</sup> V2 size standard peaks.

A sample is normalized if it is collected with a normalization size standard (specified in the primary analysis protocol [sizecalling or QC] in the assay).

**Note:** If the Sizing Quality is **X**, normalization is not applied, even if the Normalization Factor is within the normalization range. Ensure that you use the normalization size standard appropriate for your application. For more information, see "Normalization size standards provided" on page 171.

To normalize, the software:

How

applied

normalization is

- 1. Determines if the data was collected on the 3500 or 3500xL analyzer instrument.
- 2. Determines if the sample was collected with a normalization size standard definition file (normalization is enabled).
- 3. If normalization is enabled, the software calculates a Normalization Factor for the sample using multiple size standard fragments. The Normalization Factor is calculated by dividing the Normalization Target by the observed average peak height of the size standard fragments in the samples.
- 4. Compares the sample Normalization Factor to the thresholds (set in the instrument protocol).
- 5. If the calculated Normalization Factor is within the Normalization Factor range, multiplies the peak heights of the sample by the calculated Normalization Factor.

If the calculated Normalization Factor is outside the Normalization Factor range, multiplies the peak heights of the sample by the maximum or minimum Normalization Factor threshold setting (for example, if the Normalization Factor range is 0.3 to 3.0 and the calculated Normalization Factor is 5, the software applies a Normalization Factor of 3.0).

6. Indicates the normalization state of the sample in the Normalization Limit column in the Samples View.

**Normalization factor in secondary analysis** If normalization is applied in the 3500 Series Data Collection Software, the calculated Normalization factor is stored with the raw data and is applied to the raw data in the GeneMapper<sup>®</sup> *ID-X* Software v4.1 and the GeneMapper<sup>®</sup> *ID-X* Software Software v1.2 secondary analysis software. You can turn normalization off and on in the analysis method used in the GeneMapper<sup>®</sup> v4.1 and GeneMapper<sup>®</sup> *ID-X* Software v1.2 secondary analysis software. If normalization is not applied in the 3500 Series Data Collection Software (either a normalization size standard was not used, or Sizing failed  $\overbrace{}$ ), normalization cannot be applied in the secondary analysis software.

### **Review plots**

- 1. Select the samples of interest in the samples table.
- 2. Select items from the plot toolbar to manipulate the plot as needed. Place the mouse pointer over a button for the description of the button.



**IMPORTANT!** If you first view a 4-dye sample, then view a 5-dye sample, you must manually select the fifth dye. It is not automatically selected when you switch to a 5-dye sample.

Apply scaling settings to plots:
 Enter the range for Y axis and X axis, then click the Zoom buttons.

**IMPORTANT!** You must Zooming open Plot Settings each time you Zoom Y axis relative to selected peak access the View Results screen, then Zoom Y axis to 8000 click **Zoom**. Scaling settings are not automatically applied when you Zoom X axis 80 from to access this screen, or when you click hr Apply.

To apply scaling settings to all samples in the samples table, select all of the samples in the samples table to display them in the plot view, specify the scaling settings, click **Zoom**, then click **Page Up** and **Page Down** in the plot view to move through the samples.

If the plot is grayed, it indicates that the Plot Settings dialog is open. Click the 3500 task bar icon, then select Plot Settings.

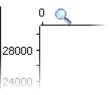
- 4. Display multiple plots as needed: in the Plot Settings Display tab, select **Checkerboard**.
- 5. Click a peak to label it (to label all peaks, see "Label peaks" on page 93).



400



- **Zoom** 1. Place the mouse pointer *above the top* of the plot or *to the left* of the plot at the start of the area you want to zoom, then click to turn the pointer to  $\bigcirc$ .
  - 2. With the still *above* the plot or to the *left* of the plot, click-drag to the end of the area you want to zoom. Do not drag the inside the plot area. Doing so changes back to a pointer and does not zoom as expected.



Change plot settings

Click (Plot Settings) in the Plot View toolbar. For information on plot settings, click (2) in the plot settings tabs.

If the *stution* is grayed, it indicates that the Plot Settings dialog is open. Click the 3500 task bar icon, then select Plot Settings.

- **Overlay samples** 1. Select samples from the Samples View to display the plots.
  - 2. Click Overlay All. When Combine Dyes is selected, the plot view displays one plot with all samples and all dyes. When Report Separate Dyes is selected, the plot view displays on plot per dye. Each dye plot contains all samples.
  - Label peaks 1. Select samples from the Samples View to display the plots.
    - 2. Click (Plot Settings) in the Plot View toolbar.



- 3. In the Plot Settings dialog box, select the Labels tab.
- 4. If you have already specified default labeling preferences, under Labelling Options:
  - a. Enable Show Peak Labels.
  - b. Click Label Peaks.
  - c. Click Apply.

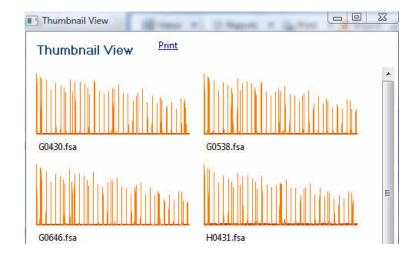
**IMPORTANT!** You must open Plot Settings each time you access the View Results screen, then click **Label Peaks**. Labelling settings are not automatically applied when you access this screen, or when you click Apply.

If you have not specified default label settings:

- a. Under Labels to Show, select the needed labels.
- b. Under Labelling Options:
  - Enable Show Peak Labels.
  - To label all peaks with the selected labels, click Label Peaks (make sure All is selected).
  - To label selected peaks, select the category from the Label Peaks list (Height, Area, Size), specify the range to label for the selected category (for example, if you select Height, specify the height range of the peaks to label), then click Label Peaks.
  - Enable **Retain Labels**.
- a. Click **Save to Preferences** to save these settings for future use. You can change preferences at any time.
- b. Click Apply.

View thumbnails Click View Thumbnails to display the traces for the samples selected in the samples view, and the dyes selected in the plot view.





### **Review sizing**

The Sizing Table View displays:

- For fragment samples All dyes
- For HID samples Size standard dye only (orange or red)

Set up the sizing	1.	Select the samples of interest in the samples table to c	lispla	y plots.	
table	2.	In the sizing table, click the Table Settings button, then specify the columns to show or hide.		Column A Table Preference Select your table display	
	3.	Filter the table as needed.	Show:	Available Columns to Show All Peaks	o Dicolau
	4.	Double-click column headers to sort columns. Multi-column sorting is supported (see "Sort" on page 97).		Show All Peaks Show Selected Peaks Show Labeled Peaks Show Selected Dye Pr	eaks
	5.	Selecting rows in the sizing table, then click Label Se	electe	ed Peaks.	
Examine the size standard plot	1.	In the Plot View toolbar, deselect all dye colors excep color (red or orange).	ot the	size standard o	lye
	2.	In the sizing table, select the size standard peaks of in	teres	t.	

3. Click Label Selected Peaks to label the size standard peaks in the Plot View.

**Note:** If labels are not displayed, click (Plot Settings) in the Plot View toolbar, then select Show Labels in the Labels tab. Click **Save to Preferences** to retain this setting.

4. Ensure that all size standard peaks are present and correctly labeled.

Overlay the sizing curve

- 1. Click (Plot Settings) in the Plot View toolbar.
- 2. Select Overlay Sizing Curve in the Display tab.

### Specify re-injections

Before the run completes, select a sample with suspect or failing flags, then click **M Re-inject**.

### View, print, and save (.pdf) sample quality reports

- 1. Select the samples of interest in the samples table.
- Click Reports to see the available reports for traces and print the reports you want.
- 3. Select the report type. Reports are displayed in the Sizing Table View at the bottom of the screen.
- 4. Modify report settings as needed.



- 5. To print the report, click 🏊 **Print**, then preview or print.
- 6. To save the report electronically (.pdf), print the report and select **CutePDF Writer** as the printer.
- 7. Close the report.



Views

🧊 Samples View

Sample Nam

Reports

View Sizing Report

View Plate Report

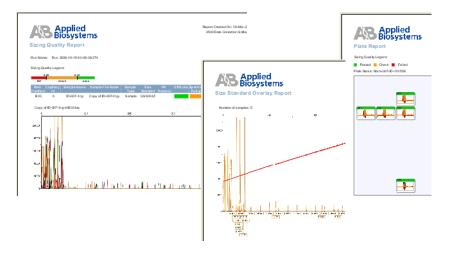
View Overlay Report

> Print

IL SIL

# • Sizing – One page per selected sample that shows the quality ranges set in the sizecalling or QC protocol, the quality values for the sample, and the electropherogram for the sample. Plot zooming is not retained in the report.

- **Overlay** One page for all selected samples that shows the size standard dyes overlaid with the size standard curves.
- **Plate** One page per plate for all selected samples that shows the well-location thumbnail traces with color-coded headers that reflect sizing quality. Plot zooming is not retained in the report.



### Export sizing results

- 1. Set up the sizing table as described above. All rows and columns displayed in the sizing table are exported.
- 2. Click Export Results.

# More features in Review Results

### **Use Rename**

**Note:** Changes to sample names are tracked only if your system includes the SAE module and auditing is enabled on your system.

- 1. In the Sample Name column, select the samples to rename, or click the Sample Name column header to select the entire column.
- 2. Click 📝 Rename.
- 3. In the Search field, enter the sample name to change.
- 4. In the Rename field, enter the new name.
- 5. Click Search, then click Rename.

### Sort

Double-click column headers to sort. Multi-column sorting is supported:

- Double-click a column header to sort the column.
- Alt+Shift-click another column header to sort another column.
- Alt+Shift-click a third column header to sort a third column.

Numbers in the column headers reflect sort order.



# Modify sequence, fragment analysis, or HID data

To edit, modify, or further analyze sequence, fragment analysis, or HID data, import the sample data files into a secondary analysis software application such as:

- Sequencing SeqScape<sup>®</sup> Software v2.7 (or later), MicroSeq<sup>®</sup> ID Analysis Software v2.2 (or later), Variant Reporter<sup>™</sup> Software (v1.1 or later), and Sequence Analysis (SeqA) Software (v5.4 or later)
- Fragment analysis GeneMapper<sup>®</sup> Software v4.1 (or later)
- **HID** GeneMapper<sup>®</sup> *ID-X* Software v1.2 (or later)

# Section 1 Calibration

## **Spatial calibration**

The 3500 Series Data Collection Software uses images collected during the spatial calibration to establish a relationship between the signal emitted by each capillary and the position where that signal falls on and is detected by the CCD camera.

### When to perform a spatial calibration

Perform a spatial calibration after you:

- Remove or replace the capillary array
- Open the detector door or move the detection cell
- Move the instrument

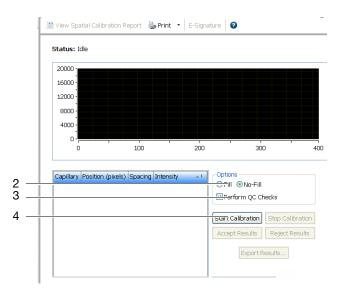
### Perform a spatial calibration

**IMPORTANT!** Do not open the instrument door during a spatial calibration run. Doing so will stop the run and require you to restart the 3500 Series Data Collection Software.

1. Access the Spatial Calibration screen: Select **Maintenance**, then select **Spatial Calibration** in the navigation pane.

**Note:** The screen does not display results unless you have previously performed a spatial calibration.





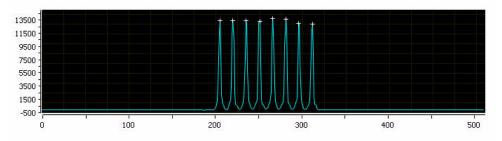
2. Select **No Fill**, or select **Fill** to fill the array with polymer before starting the calibration.

(Optional) Select **Perform QC Checks** if you want the system to check each capillary against the specified range for spacing and intensity. During the calibration, the software calculates:

Attribute	Calculation	Threshold
Average peak height	sum of all peak heights number of peaks	<ul><li>8-cap: 6400 RFU</li><li>24-cap: 3000 RFU</li></ul>
Uniformity (peak height similarity)	standard deviation average peak height	0.2
Capillary spacing	max spacing - min spacing	2 pixels

#### 3. Click Start Calibration.

The display updates as the run progresses.



If the average of any of the QC values exceeds the threshold, a Spatial QC Check error message is displayed.

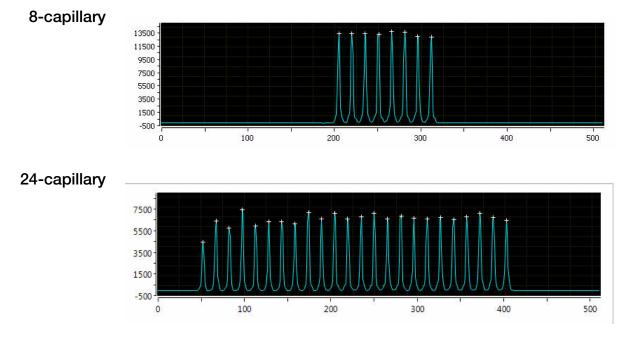
### Evaluate the spatial calibration profile

When the run is complete:

- 1. Evaluate the spatial calibration profile to ensure that you see:
  - One sharp peak for each capillary. Small shoulders are acceptable.
  - One marker (+) at the apex of every peak. No off-apex markers.
  - An even peak profile (all peaks about the same height).
- 2. If the results meet the criteria above, click Accept Results.

If the results do not meet the criteria above, click **Reject Results**, then go to "Spatial calibration troubleshooting" on page 300.

### Example spatial profiles



### Export spatial calibration results

To export spatial calibration results:

- 1. Click Export.
- 2. Enter an export file name.



4. Click Save.

Options ○ Fill				
Start Calibration	Stop Calibration			
Accept Results	Reject Results			
Export Results				

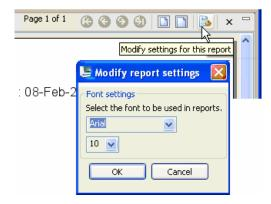
The export file contains the following results:

- Capillary Number
   Spacing
- Position (pixels)
   Intensity

### View and print a spatial calibration report

**Note:** Spatial and spectral calibration reports include the date on which a capillary array is installed for the first time on the instrument. Install standard reports use the most recent install date if a capillary array was removed and re-installed on the instrument.

- 1. Click 📰 View Spatial Calibration Report.
- 2. In the Report screen, click toolbar options to manipulate the report as needed. Place the mouse pointer over an item for a description of the item.



3. To print the report, click 🏊 **Print**.



### Save historical calibration reports (.pdf) for record keeping

**IMPORTANT!** After performing a calibration, save the calibration report electronically for record keeping. The software does not save historical calibration results. Only the most recent spatial calibration is maintained in the software.

- 1. Click 🔛 View Spatial Calibration Report.
- 2. Click 🌺 Print.

4. Close the report.

- 3. In the Printer dialog box, select CutePDF Writer as the printer.
- 4. Specify a name and location for the report.

## **Spectral calibration**

A spectral calibration creates a de-convolution matrix that compensates for dye overlap (reduces raw data from the instrument) in the 4-dye, 5-dye, 6-dye, or AnyDye data stored in each sample file.

### When to perform a spectral calibration

Perform a spectral calibration for each dye set/polymer type combination you will use:

- Sequencing dye set/polymer type
- Fragment dye set/polymer type
- HID dye set/polymer type

Perform a spectral calibration when you:

- Use a dye set that you have not previously calibrated
- Change the capillary array
- Change the polymer type
- Have a service engineer perform an optical service procedure, such as realigning or replacing the laser or CCD camera or mirrors on the instrument
- See a decrease in spectral separation (pull-up/pull-down in peaks) in the raw or analyzed data

**Note:** If you are using the v3.1 sequencing standard or v1.1 sequencing standard and want to run a performance check and a spectral calibration, you can skip this process, and run the Sequencing Install Standard performance check. If you select Keep Spectral Calibration Data in the Performance Check, the software runs a spectral calibration for dye set E or Z during a sequencing check and allows you to save the spectral calibration data. For information, see "Run the sequencing install standard performance check" on page 119.

# Estimated run times

Standard	Polymer Type	Run Time (min)
Matrix standard	Any	≤30
Sequencing standard	POP-7 <sup>™</sup> polymer	≤40
	POP-6 <sup>™</sup> polymer	≤135

### Prepare for the spectral calibration

Prepare the<br/>instrument1. If you have not already done so, perform a spatial calibration (see "Spatial<br/>calibration" on page 99).

- 2. In the Dashboard, check consumable status (page 29). Ensure that:
  - Consumables are not expired
  - Adequate injections remain for consumables
- 3. Ensure that the buffer levels are at the fill lines ("Check buffer fill levels" on page 31).
- 4. Set the oven temperature, then click **Start Pre-heat**:
  - **60** °C POP-7<sup>TM</sup> polymer
  - **50** °C POP-6<sup>TM</sup> polymer

Pre-heat the oven and detection cell while you prepare for a run (detection cell temperature is set by the software). Preheating helps mitigate subtle first-run migration rate effects. The pre-heat function automatically turns off after 2 hours.

Applied Biosystems recommends that you pre-heat the oven for at least 30 minutes before you start a run if the instrument is cold.

5. Check the pump assembly for bubbles and run the Remove Bubble wizard if needed (see page 251).

Prepare the standard calibration plate **IMPORTANT!** Do not use warped or damaged plates.

1. Prepare the calibration standard as described in the standard product insert. See Table 28 on page 259 and Table 29 on page 260 for standard part numbers.

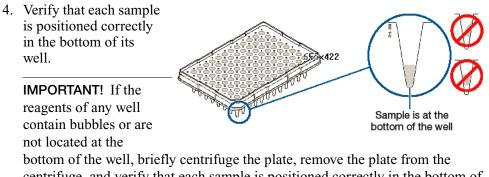
Dye set	Standard							
Е	BigDye® Terminator (BDT) v1.1 Sequencing Standard							
	BigDye® Terminator (BDT) v1.1 Matrix Standard							
Z	BigDye® Terminator (BDT) v3.1 Sequencing Standard							
	BigDye® Terminator (BDT) v3.1 Matrix Standard							
F	DS-32 Matrix Standard							
E5	DS-02 Matrix Standard							
G5	DS-33 Matrix Standard							

2. Load the standards in injection position 1 in the spectral calibration plate:

**IMPORTANT!** You do not create a plate for the calibration. The software uses predetermined positions for the calibration. You cannot specify standard location on the plate. If you do not place calibration standards in the positions indicated, the calibration will fail.

8-capillary	A1 through H1
96-well plate	
24-capillary	A1 through H1, A2 through H2, and
1 2	A3 through H3
96-well plate	
	4
24-capillary	Columns 1, 3, vozsraumian monos
1 2	and 5 in rows
384-well plate	A, C, E, G, I, K,
Note: 384-well	M, O
plates are not	
supported on 8-capillary	
instruments.	

- **96** Supports 96-well standard reaction plate. 8-strip standard tubes are also supported with appropriate retainers.
- **96-Fast Tube** Supports 96-well Fast reaction plate. 8-strip fast tubes are also supported with appropriate retainers
- 3. Briefly centrifuge the plate containing the standards.



centrifuge, and verify that each sample is positioned correctly in the bottom of its well.

5. Store the plate on ice until you prepare the plate assembly and load the plate in the instrument.

#### Prepare the plate **IMPORTANT!** Prepare the plate assembly on a clean, level surface. Do not heat plates that are sealed with septa. assembly

- 1. Align the holes in the septa strip with the wells of the plate, then firmly press downward onto the plate.
- 2. Place the sample plate into the plate base.

Plate retainer Plate with septa strip **IMPORTANT!** Make sure to use the correct plate base for Plate base standard plates versus 8-tube strips and fast plates. Using the wrong plate base may affect

- 3. Snap the plate retainer (cover) onto the plate, septa, and plate base.
- 4. Verify that the holes of the plate retainer and the septa strip are aligned. If not aligned, re-assemble and then assemble the plate assembly.

**IMPORTANT!** The array tips will be damaged if the plate retainer and septa strip holes do not align correctly.

#### Load the plate in 1. Place the plate in the autosampler with the labels the instrument facing you (or the instrument door) and the notched corner of the plate in the notched corner of the autosampler.

performance.

- 2. Close the instrument door to re-initialize the instrument.

### Perform a spectral calibration

**IMPORTANT!** Do not change electronic signature settings during a spectral calibration.

**IMPORTANT!** If you change polymer type, spectral calibrations for the original polymer type are not retained.

 Access the Spectral Calibration screen: Select Maintenance, then select Spectral Calibration in the navigation pane.

**Note:** The screen does not display results until you perform a spectral calibration. To view previous calibration data, click **History View**.

Library	Maintenance	Tools	1	Manage	
	3				
Calibra	te		1		
		Spatia	8		
		Spectra	1/		
Devterme	ince check				

Calibration Settings									0	Current In: Polymer Ty	strument Consuma pe: POP4	bles Capillary I	.ength
Number of Wells: (	96 🔿	96-Fast	tTube	0 384		Cher	nistry S	tandard:	<b></b>		Start Run	]	
Plate Position: 🤇	A OB						[	Dye Set:	~				0%
Allow Borrowing										Status: F	Ready		
▼ Capillary Run Da	ta												
Capillary	1	2	3	4	5	6	7	8					
Run 1													
Run 2													
Run 3													
Overall													
Passed	Faile	ed		Borrow	ed		Not Cal	ibrated					
		Qua	lity Va	lue:			Con	dition #:	Status:		Message:		
▼ Intensity vs Sca	n Numbe	r											
Raw Da	a	*						<b>~</b>	] 🗹 🗹 🗹 🗹			33(+)	3 <b>0</b>
0	4000	)		8000		1200	0	1600	0 20000	24000	28000	32000	1.101.00
1									· · · · ·		· · · · ·	1	
40000 -													
20000 -													
-													
0													
								Intensity v	s Scan Number				

2. Select the number of wells in the spectral calibration plate and specify the plate location in the instrument.

**Note:** You do not create a plate for the calibration. The software uses predetermined positions for the calibration. You cannot specify standard location on the plate.

3. Select the chemistry standard and the dye set that you are running the calibration for.

**Note:** If the dye set list is empty, ensure that your instrument is configured with a compatible polymer type and capillary length for the selected chemistry standard.

**IMPORTANT!** To calibrate a custom dye set using AnyDye, first create the dye set (see "Create a new dye set" on page 168), then select the name of the custom dye set from the Dye Set list. The AnyDye selection in the Dye Set list contains default settings. It does not correspond to custom dye sets created with the AnyDye dye set template.

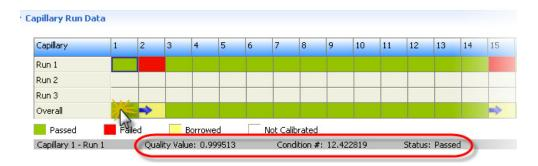
- 4. (Optional) Select **Allow Borrowing**. Selecting this option instructs the software to automatically replace information from failed capillaries with information from an adjacent passing capillary with the highest Quality value. For more information, see "What you see during a spectral calibration" on page 112.
- 5. Click Start Run. The following occurs:
  - The system sets up three injections (see "What you see during a spectral calibration" on page 112 for information on the number of injections performed).
  - The Capillary Run Data display updates after each injection is complete.
  - The status bar updates during Run 1.

**IMPORTANT!** The status bar does not update during Run 2 or Run 3.

• Passing and failing capillaries are shown in green and red respectively. Borrowed capillaries are shown in yellow with an arrow indicating the adjacent capillary from which results were borrowed.

To display the result for each capillary (spectral data, Quality Value, and Condition Number) below the run results table, click a capillary in the table.

**Note:** The results displayed when you click a borrowed capillary are the passing results borrowed from the adjacent capillary. To determine the reason that a capillary fails, view the spectral calibration report. See "View and print a spectral calibration report" on page 116.



For all spectral calibration injections (even capillaries that are green in the Overall row), evaluate the data as described in the next section.

Spectral Quality Values and Condition Numbers

#### Spectral Quality Value

A spectral Quality Value reflects the confidence that the individual dye emission signals can be separated from the overall measured fluorescence signal. It is a measure of the consistency between the final matrix and the data from which it was computed. A Quality Value of 1.0 indicates high consistency, providing an ideal matrix with no detected pull-up/pull-down peaks.

In rare cases, a high Quality Value can be computed for a poor matrix. This can happen if the matrix standard contains artifacts, leading to the creation of one or more extra peaks. The extra peak(s) causes the true dye peak to be missed by the algorithm, and can lead to a higher Quality Value than would be computed with the correct peak. Therefore, it is important to visually inspect the spectral calibration profile for each capillary (see "Evaluate the spectral calibration data" on page 110).

#### **Condition Number**

A Condition Number indicates the amount of overlap between the dye peaks in the fluorescence emission spectra of the dyes in the dye set.

If there is no overlap in a dye set, the Condition Number is 1.0 (ideal conditions), the lowest possible value. The condition number increases with increasing peak overlap.

The ranges that the software uses to determine if a capillary passes or fails are:

Dye Set	Quality Value Minimum	Condition Number Maximum
AnyDye	0.8 (default)	20.0 (default)
E	0.95	5.5
E5	0.95	6.0
F	0.95	8.5
G5	0.95	13.5
J6	0.95	8.0
Z	0.95	5.5

### Evaluate the spectral calibration data

**IMPORTANT!** Do not accept a spectral calibration until you examine the data for all capillaries.

When a spectral calibration completes successfully, the Overall row displays green, red, or yellow results.

For each capillary:

- 1. Click a capillary to display the spectral and raw data for a capillary.
- 2. Check that the data meet the following criteria:

Attribute	Acceptance Criteria	Example
Order of the peaks in the spectral profile from left to right	4-dye: blue-green-yellow-red	Elue Green Yellow Red
	• 5-dye: blue-green-yellow-red-orange	Blue Groen Yellow Red Orange
Order of the peaks in the raw data profile from left to right	<ul> <li>Sequencing (matrix standard only):</li> <li>4-dye: red-yellow-blue-green</li> </ul>	Ped Yillow Blue Green
	<ul> <li>Fragment analysis/HID:</li> <li>4-dye: red-yellow-green-blue</li> <li>5-dye: orange-red-yellow-green-blue</li> </ul>	Oringe Bed Yelkow Green Blue
Extraneous peaks in the raw data profile	None <b>Note:</b> The E5 profile may include extraneous peaks outside the matrix peak region which can be ignored.	

Attribute	Acceptance Criteria	Example
Peak morphology in the spectral profile	<ul> <li>No gross overlaps, dips, or other irregularities</li> <li>Peaks separate and distinct</li> </ul>	
	<b>Note:</b> The profiles of G5 (shown to the right), F, and J6 may not be as smooth as the profiles for other dye sets (shown above) due to the effect of variable binning (a feature that reduces signal variation between dyes of different fluorescent efficiencies).	

- 3. As needed, zoom on the spectral profile traces to determine if the data meet the criteria:
  - a. Place the pointer *above the top* of the plot or *to the left* of the plot at the start of the area you want to zoom, then click to turn the pointer to  $\bigcirc$ .



b. With the still *above* the plot or to the *left* of the plot, click-drag to the end of the area you want to zoom. Do not drag the sinside the plot area. Doing so changes back to a pointer and does not zoom as expected.

You can also click zoom and fit buttons to zoom. 🐑 腔

- 4. If the data for all capillaries meet the criteria above, click Accept Results.
- 5. If any capillary data does not meeting the criteria\_above, click **Reject Results**, then go to "Spectral calibration troubleshooting" on page 301.

### What you see during a spectral calibration

A spectral calibration automatically sets up three injections. The number of injections performed depends on:

- The number of capillaries that pass or fail during an injection
- Whether you select the Allow Borrowing option

**Note:** The first time you perform a spectral calibration (for each dye set) after installing a new capillary array, you may notice pull-down peaks (or mirror image peaks). These pull-down peaks will eventually correct themselves once the run completes.

Capillary information sharing

Spectral

calibration with

A spectral calibration can share capillary information:

- **Between injections** If a capillary in an injection does not meet the spectral Quality Value and Condition Number limits shown on page 109, the software automatically uses the information from that capillary in a different injection.
- Within an injection If a capillary in an injection does not meet the spectral Quality Value and Condition Number limits shown on page 109 and the Allow Borrowing option is selected, the software can also use the information from a capillary to the left or the right of that capillary, if the values are higher than those for that capillary in a different injection.

When Borrowing is *disabled*, all capillaries must pass (meet the spectral Quality Value and Condition Number limits) for the calibration to pass.

Jan		value and C	ondition realiser minus) for the carloration to pass.
	Borrowing disabled	Injection 1	The software evaluates the Quality Value and Condition Number of all capillaries.
			<ul> <li>If all capillaries pass, the calibration is complete, and injections 2 and 3 are not performed.</li> </ul>
			If any capillaries fail, injection 2 is performed.
		Injection 2	• The software evaluates the Quality Value for each capillary across injections 1 and 2 and uses the information from the capillary with the highest Quality Value.
			• If all capillaries now pass, the calibration is complete and injection 3 is not performed.
			<ul> <li>If the same capillary fails in both injection 1 and 2, injection 3 is performed.</li> </ul>
		Injection 3	• The software evaluates the Quality Value for each capillary across injections 1, 2, and 3 and the information from the capillary with the highest Quality Value.
			<ul> <li>If all capillaries now pass, the calibration passes.</li> </ul>
			• If the same capillary fails in injection 1, 2, or 3, the calibration fails.

### Spectral calibration with Borrowing enabled

Allow Borrowing

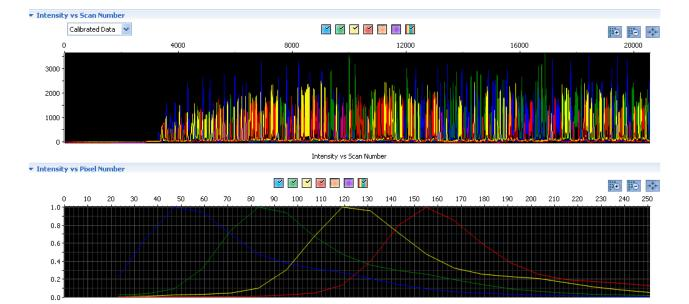
When Borrowing is *enabled*, all capillaries have to pass (meet the spectral Quality Value and Condition Number limits) within the borrowing limits:

- 8-capillary instruments One adjacent-capillary borrowing event allowed
- 24-capillary instruments Up to three adjacent-capillary borrowing events allowed (the number of allowed borrowing events can be decreased in Preferences).

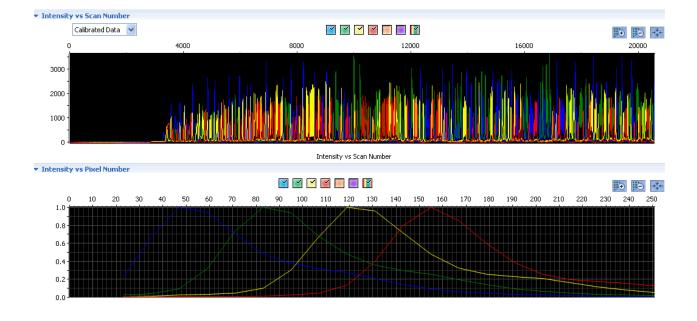
Injection 1	<ul> <li>The software evaluates the Quality Value and Condition Number of all capillaries.</li> </ul>
	<ul> <li>If all capillaries pass, the calibration is complete, and injections 2 and 3 are not performed.</li> </ul>
	• If any capillaries fail, the software borrows from an adjacent capillary.
	• If, after borrowing, >1 or > 3 capillaries fail, injection 2 is performed.
Injection 2	• The software evaluates the quality values between adjacent capillaries in injection 2 and for each capillary across injections 1 and 2 and the information with the highest Quality Value for each capillary.
	<ul> <li>If all capillaries pass, the calibration is complete and injection 3 is not performed.</li> </ul>
	<ul> <li>If, after borrowing, &gt;1 or &gt; 3 capillaries from injection 1 or 2 do not pass, injection 3 is performed.</li> </ul>
Injection 3	• The software evaluates the quality values between adjacent capillaries in injection 3 and for each capillary across injections 1, 2, and 3, then the information with the highest Quality Value for each capillary.
	<ul> <li>If all capillaries now pass, the calibration passes.</li> </ul>
	<ul> <li>If after borrowing, &gt;1 or &gt; 3 capillaries from injection 1, 2, or 3 do not pass, the calibration fails.</li> </ul>
	1

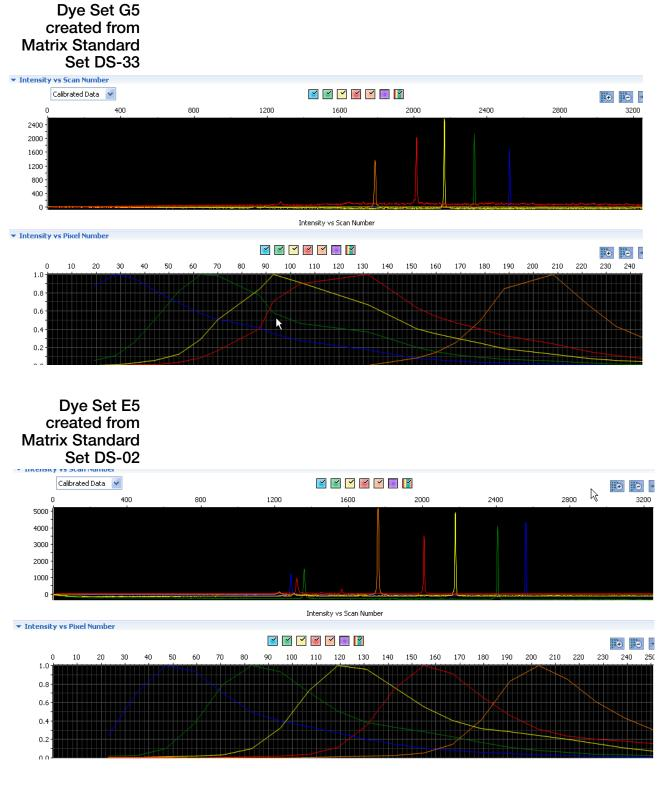
### Example spectral calibration data

Dye Set E created from Sequencing Standard



### Dye Set Z created from Sequencing Standard





### Export spectral calibration results

To export spectral calibration results:

1. Click 🛃 Export Spectral Calibration Results.

- 2. Specify an export file name and location.
- 3. Click Save.

The export file contains the following results:

- Capillary Number
- Condition Number
- Scan Number

- Quality Value
- Peak Height
- Reason For Failure
- Borrowed From Capillary Run From Injection

### View and print a spectral calibration report

**Note:** Spatial and spectral calibration reports include the date on which a capillary array is installed for the first time on the instrument. Install standard reports use the most recent install date if a capillary array was removed and re-installed on the instrument.

- 1. Click 📰 View Spectral Calibration Report.
- 2. In the Report screen, click toolbar options to manipulate the report as needed. Place the mouse pointer over an item for a description of the item.

Page 1 of 1	C C O O O D	
	Modify report settings 🛛	
:08-Feb-2	<ul> <li>Font settings</li> <li>Select the font to be used in reports.</li> <li>Aria</li> <li>IO</li> </ul>	
	OK Cancel	
3. To print	the report, click Ъ Prin	t.
4. Close th	e report. Page 1 of 1	6666

### Save historical calibration reports (.pdf) for record keeping

**IMPORTANT!** After performing a calibration, save the calibration report electronically for record keeping. The software does not save historical calibration results. Only the most recent spectral calibration for each dye set is maintained in the software.

- 1. Click 📄 View Spectral Calibration Report.
- 2. Click 🏊 Print.
- 3. In the Printer dialog box, select **CutePDF Writer** as the printer.
- 4. Specify a name and location for the report.

### View the spectral calibration history

Select History View, then select a dye set to view the associated calibration history.

Calibration	n Information 🥖 🚬		
Dye Set	Chemistry Standard	Calibration Date 🗸 🗸	Capillary Arro
F	Matrix Standard	20-Jan-2009 11:38:52 AM	80K2450
G5 💥	Matrix Standard	13-Jan-2009 10:11:36 PM	80K2450

# Section 2 Performance check

The Performance check allows you to periodically self-check the instrument system using Applied Biosystems standard.

### Run the sequencing install standard performance check

**When to perform** When your instrument is installed, the service engineer runs a sequencing install standard performance check.

Applied Biosystems recommends that you run the sequencing install standard performance check monthly to verify that the instrument meets read length specifications.

The Sequencing Install Performance check has an option to include and save the spectral calibration. If you select this option and you accept the sequencing install standard results, you do not need to run the spectral calibration (described in "Spectral calibration" on page 103) for E and Z dye sets. You still need to run spectral calibrations for other dyes sets.

The performance check is application-specific. If you will run general sequencing applications with POP-7<sup>TM</sup> polymer and MicroSeq<sup>®</sup> ID applications with POP-6<sup>TM</sup> polymer, install the appropriate polymer and perform separate performance checks.

Estimated run	• General sequencing – 45 minutes
times	• MicroSeq <sup>®</sup> ID $- 2$ hours

### Prepare for the sequencing install standard performance check

Prepare the	1. In the Dashboard, check consumable status ("Check consumable status" on
instrument	page 29). Ensure that:

- Consumables are not expired
- Adequate injections remain for consumables
- 2. Ensure that the buffer levels are at the fill lines ("Check buffer fill levels" on page 31).
- 3. Set the oven temperature, then click Start Pre-heat:
  - **60** °C General sequencing POP-7<sup>TM</sup> polymer
  - **50** °C MicroSeq<sup>®</sup> ID POP-6<sup>TM</sup> polymer

Pre-heat the oven and detection cell while you prepare for a run (detection cell temperature is set by the software). Preheating helps mitigate subtle first-run migration rate effects. The pre-heat function automatically turns off after 2 hours.

Applied Biosystems recommends that you pre-heat the oven for at least 30 minutes before you start a run if the instrument is cold.

4. Check the pump assembly for bubbles and run the Remove Bubble wizard if needed (see page 251).

Prepare the installation standard plate **IMPORTANT!** Do not use warped or damaged plates.

- $\bigcirc$
- 1. Prepare the sequencing install standard as described in the product insert. See Table 28 on page 259 for standard part numbers.

Application	Standard
General sequencing (POP-7 <sup>™</sup> polymer, 50-cm capillary)	BigDye <sup>®</sup> Terminator (BDT) v3.1 Standard
MicroSeq <sup>®</sup> ID applications (POP-6 <sup>™</sup> polymer, 50-cm capillary)	BigDye <sup>®</sup> Terminator (BDT) v1.1 Standard

2. Load the standards in injection position 1 in the spectral calibration plate:

**IMPORTANT!** You do not create a plate for the performance check. The software uses predetermined positions for the performance check run. You cannot specify standard location on the plate. If you do not place standards in the positions indicated, the calibration will fail.

8-capillary	A1 through H1
96-well plate	
24-capillary	A1 through H1, A2 through H2, and
96-well plate	A3 through H3
·	
	<b>ω</b>
24 appillant	Columna 1.2
24-capillary	Columns 1, 3, and 5 in rows
384-well plate	A, C, E, G, I, K,
Note: 384-well	M, O
plates are not	
supported on 8-capillary	
instruments.	0

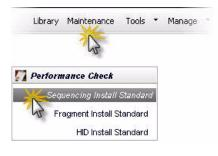
- **96** Supports 96-well standard reaction plate. 8-strip standard tubes are also supported with appropriate retainers.
- **96-Fast Tube** Supports 96-well Fast reaction plate. 8-strip fast tubes are also supported with appropriate retainers.
- 3. Briefly centrifuge the plate containing the standards.

	4. Verify that each sample is positioned correctly in the bottom of its well. IMPORTANT! If the reagents of any well contain bubbles or are not located at the bottom of the well, briefly centrifuge the plate, remove the plate from the centrifuge, and verify that each sample is positioned correctly in the bottom of its well.
	5. Store the plate on ice until you prepare the plate assembly and load the plate in the instrument.
Prepare the plate assembly	<b>IMPORTANT!</b> Prepare the plate assembly on a clean, level surface. Do not heat plates that are sealed with septa.
	1. Align the holes in the septa strip with the wells of the plate, then firmly press downward onto the plate.
	2. Place the sample plate into the plate base. Plate with septa strip
	IMPORTANT! Make sure to use the correct plate base for standard plates versus 8-tube strips and fast plates. Using the wrong plate base may affect performance.
	3. Snap the plate retainer (cover) onto the plate, septa, and plate base.
	4. Verify that the holes of the plate retainer and the septa strip are aligned. If not aligned, re-assemble and then assemble the plate assembly.
	<b>IMPORTANT!</b> The array tips will be damaged if the plate retainer and septa strip holes do not align correctly.
Load the plate in the instrument	<ol> <li>Place the plate in the autosampler with the labels facing you (or the instrument door) and the notched corner of the plate in the notched corner of the autosampler.</li> </ol>
	2. Close the instrument door to re-initialize the

2. Close the instrument door to re-initialize the instrument.

### Run the sequencing install standard performance check

- Access the sequencing install standard screen (Figure 9 on page 122): Select Maintenance, then select Sequencing Install Standard in the navigation pane.
- 2. Select the chemistry type: General Sequencing or MicroSeq<sup>®</sup> ID.



3. Select the plate type and plate position in the instrument.

**Note:** You do not create a plate for the performance check. The software uses predetermined positions for the run. You cannot specify standard location on the plate.

CRL Pass/Fail Threshold:       CCRL Pass/Fail Threshold:       Current Instrument Consumables         Creating v       Read Length Start:       Polymer Type:       POlymer Typ	Calbration Settings			Crosing	Catting											
CRL Pass/Fail Threshold: Sool S   Polymer Type: Polymer Type:	Calloration Develops						_		Curren	t Instrum	ent Con	simable				
Image: solution of Wells: 0.96 0.96-FastTube 0.384       Read Length Start: 400 0.0000       Read Length Start: 400 0.0000         Plate Position: 0.A 0.8       Read Length End: 539 0.0000       Status: Ready         Image: solution Data       Read Length End: 539 0.0000       Status: Ready         Image: solution Data       Read Length End: 539 0.0000       Status: Ready         Image: solution Data       Image: solution Data       Image: solution Data         Image: solution Data       Image: solution Data       Image: solution Data         Image: solution Data       Image: solution Data       Image: solution Data         Image: solution Data       Image: solution Data       Image: solution Data         Image: solution Data       Image: solution Data       Image: solution Data         Image: solution Data       Image: solution Data       Image: solution Data         Image: solution Data       Image: solution Data       Image: solution Data         Image: solution Read Length (CRL)       Image: solution Data       Image: solution Data         Image: solution Read Length (CRL)       Image: solution Data       Image: solution Data         Image: solution Read Length       Image: solution Data       Image: solution Data         Image: solution Read Length (CRL)       Image: solution Read Length       Image: solution Data         Image:				CRL Pa	ss/Fail T	hreshold	500							y Length	: 36cm	n
Image: solution of Wells: 0.96 0.96-FastTube 0.384       Read Length Start: 400 0.0000       Read Length Start: 400 0.0000         Plate Position: 0.A 0.8       Read Length End: 539 0.0000       Status: Ready         Image: solution Data       Read Length End: 539 0.0000       Status: Ready         Image: solution Data       Read Length End: 539 0.0000       Status: Ready         Image: solution Data       Image: solution Data       Image: solution Data         Image: solution Data       Image: solution Data       Image: solution Data         Image: solution Data       Image: solution Data       Image: solution Data         Image: solution Data       Image: solution Data       Image: solution Data         Image: solution Data       Image: solution Data       Image: solution Data         Image: solution Data       Image: solution Data       Image: solution Data         Image: solution Data       Image: solution Data       Image: solution Data         Image: solution Read Length (CRL)       Image: solution Data       Image: solution Data         Image: solution Read Length (CRL)       Image: solution Data       Image: solution Data         Image: solution Read Length       Image: solution Data       Image: solution Data         Image: solution Read Length (CRL)       Image: solution Read Length       Image: solution Data         Image:	General Sequer	cing 🗸										_				
Plate Position: <ul> <li></li></ul>	tumber of Wells: 0 96 0 96-Fa	stTube 038	4	R	ead Leng	gth Start	40 🗘					Sta	art Run	J		
	Plate Position:    A							_					_			
Capillary Run Data      Contiguous Read Length (CRL)      CRL Pass/Fail      Comparison with Ref Sequence      CRL Basepair Accuracy      Basepair Accuracy      Read Length      Passed      Failed      Borrowed      Not Calibrated      Condition #:      Status:      Message:      Intensity vs Scan Number      Raw Data      Yo Scan Number      Raw Data      Yo Scan Number      Read Length      Yo	Spectral Calibration Data			1	Read Ler	ngth End	539		Statu	is: Read	У					
Capillary       1       2       3       4       5       6       7       8       9       10       11       12       13       14       15         Spectral Calibration Run																
Spectral Calbration Run Contiguous Read Length (CRL) CRL Pass/Fail Comparison with Ref Sequence CRL Basepair Accuracy Ba	▼ Capillary Run Data												_1			
Spectral Calbration Run Contiguous Read Length (CRL) CRL Pass/Fail Comparison with Ref Sequence CRL Basepair Accuracy Ba			_	_	_		-	_	_	_	_	_		_	_	_
Contiguous Read Length (CRL) CRL Pass/Fail Comparison with Ref Sequence CRL Basepair Accuracy Basepair A		1	2	3	4	5	6	7	8	9	10	11	12	13	14	15
CRL Pass/Fail																
Comparison with Ref Sequence     Image: Comparison with Ref Sequence     Image: Comparison with Ref Sequence       CRL Basepair Accuracy     Image: Comparison with Ref Sequence     Image: Comparison with Ref Sequence       CRL Basepair Accuracy     Image: Comparison with Ref Sequence     Image: Comparison with Ref Sequence       Basepair Accuracy     Image: Comparison with Ref Sequence     Image: Comparison with Ref Sequence       Passed     Faled     Borrowed       Quality Value:     Condition #:     Status:       Message:     Image: Condition #:     Status:       Intensity vs Scan Number     Image: Condition #:     Image: Comparison #:       Raw Data     Image: Comparison #:     Image: Comparison #:				_	_			-			_		_			-
CRL Basepair Accuracy Basepair Accuracy Basepair Accuracy Read Length  Passed Failed Borrowed Quality Value: Condition #: Status: Message:  Intensity vs Scan Number  Raw Data										_	_					
Basepair Accuracy     Read Length       Passed     Failed       Passed     Failed       Borrowed     Not Calbrated       Condition #:     Status:       Message:		ce			_											
Read Length     Image: Condition #1     Status:     Message:       Passed     Faled Qualky Value:     Condition #1     Status:     Message:       Intensity vs Scan Number     Image: Condition #2     Image: Condition #2     Image: Condition #2       Raw Data     Image: Condition #2     Image: Condition #2     Image: Condition #2		_			_		_			_						
Passed ■ Faled Borrowed Not Calibrated     Quality Value: Condition #: Status: Message:      Intensity vs Scan Number     Raw Data      Passed ■ Passe																
Passed     Faled     Borrowed     Not Calibrated       Quality Value:     Condition #:     Status:     Message:       ✓ Intensity vs Scan Number     Raw Data     ✓     ✓     ✓	Read Length															
Quality Value:     Condition #:     Status:     Message:       Intensity vs Scan Number     Raw Data     Image: Condition #:     Status:     Message: Condition #:																
▼ Intensity vs Scan Number Raw Data ▼			d	N					0.0	ha a ma						_
Raw Data 👻		alky value;			Con	aloon #:	_		268	cus:			P	essage:		
								_								
0 4000 8000 12000 16000 20000 24000 28000 32000	Raw Data 💌						1 🗹 🗅	1 🗹	<b>I</b>							۲
	0 4000	8000		12	000		16000		20000		2400	0	28	000	3	2000
	Ì															
	0.															
0-																
						Inten	sity vs Sc	an Nur	nber							
Intensity vs Scan Number	Sequence Comparison to Sam	ple (Capillary	(1)													

Figure 9 Sequencing Install Standard screen

4. (Optional) If you have not already run a spectral calibration, select **Keep Spectral Calibration Data** to save the sequencing install standard run (if it passes) as a spectral calibration. General Sequencing with BDTv3.1 Install standard and POP-7<sup>™</sup> polymer generates a Z dye set spectral calibration.

**Note:** he spectral calibration record will only be saved if Keep Spectral Calibration Data option is checked on the screen. If you decide to uncheck the option, create a separate spectral calibration from the Maintenance menu.

- MicroSeq<sup>®</sup> with BDTv1.1 Install Standard and POP-6<sup>™</sup> polymer generates an E dye set spectral calibration.
- 5. Click Start Run.

**IMPORTANT!** Do not accept a sequencing installation standard run until you examine the data.

### What you see during a run

The system performs one run, then evaluates:

- Spectral data, if you specified to keep spectral data
- Sequence data

The Capillary Run Data display (Figure 10 on page 124) updates after the run is complete:

• The spectral calibration status is displayed in the first row of the run results table. Passing and failing capillaries in the performance run are shown in green and red respectively for the CRL criteria. Borrowed capillaries (spectral only) are shown in yellow with an arrow indicating the adjacent capillary from which results were borrowed. The spectral result for each capillary is displayed below the run results table.

**Note:** Clicking a borrowed capillary displays borrowed, not failed, data. For information on why a capillary failed, look in the Sequencing Install Standard Detail Report.

- The sequencing install standard status is displayed in the third row of the run results table (CRL Pass/Fail).
- The Quality Value and Condition Number for each capillary is displayed below the table.

Capillary	1	2	3	4	5	6	7	8	9	10	Median	SD
Spectral Calibration Run			-									
Contiguous Read Length (CRL)	691.0	712.0	0.0	711.0	705.0	703.0	708.0	704.0	704.0	710.	708.5	144.98
CRL Pass/Fail												
Comparison with Ref Sequence												
CRL Basepair Accuracy	100.0	99.9		99.9	99.9	99.9	99.9	99.9	99.9	99.9	99.9	
Basepair Accuracy	100.0	100.0		100.0	100.0	100.0	100.0	100.0	100.0	100.	100.0	
Read Length	500.0	500.0		500.0	500.0	500.0	500.0	500.0	500.0	500.	500.0	102.47

Figure 10 Sequencing install standard – capillary run data

# How the software determines passing and failing capillaries for the spectral calibration

The software evaluates the Quality Value and Condition Number for each capillary (for more information, see "Spectral Quality Values and Condition Numbers" on page 109).

Borrowing is automatically enabled: 1 borrowing event is allowed for 8-capillary instruments, up to 3 borrowing events for 24-capillary instruments. For more information, see "Capillary information sharing" on page 112. The number of borrowing events can be decreased – see "User preferences" on page 34.

Dye Set	Quality Value Minimum	Condition Number Maximum
E	0.95	5.5
Z	0.95	5.5

Thresholds used by the software for pass/fail are:

# How the software determines passing and failing capillaries for the sequencing performance check

The software calculates the Contiguous Read Length for each capillary. Capillaries that are below the threshold fail. The remaining results that the software displays are for information only.

	Description
Contiguous Read Length (CRL)	The longest uninterrupted segment of bases with an average Quality Value (QV) $\ge$ 20.
	In addition to evaluating the QV of a base call, the software considers the QV of adjacent bases within a $\pm$ 20-bp moving average to determine a contiguous read length based on quality values: the software starts from the 5' end and calculates the average QV across a moving window size of 20, sliding 1 bp at a time, to the 3' end. The resulting longest contiguous segment is determined as the CRL.
CRL Pass/Fail	• General sequencing – Capillaries with a CRL ≤ 500 bp fail.
	• <b>MicroSeq<sup>®</sup> ID</b> – Capillaries with a CRL $\leq$ 600 bp fail.
For information only sequencing install star	- Based on alignment of the base-called sample sequence with the known reference of the ndard
CRL Basepair	CRL accuracy is determined by base-pair comparison between the base-called sample and
Accuracy	the known reference sequence for the install standard within the contiguous read length region calculated (as described in the CRL definition above).
	the known reference sequence for the install standard within the contiguous read length
Accuracy	the known reference sequence for the install standard within the contiguous read length region calculated (as described in the CRL definition above).
Accuracy	<ul> <li>the known reference sequence for the install standard within the contiguous read length region calculated (as described in the CRL definition above).</li> <li>The length of read (in bases) at which base calling accuracy is ≥98.5%.</li> <li>The read length value for this information is derived from basecall-accuracy, not from quality</li> </ul>

### Evaluate sequencing install standard data

When a sequencing install standard run completes successfully, the CRL Pass/Fail row displays green or red results.

For each capillary:

- 1. Click a capillary to display the spectral and raw data profiles for a capillary.
- 2. Check that the data meet the following criteria:

Attribute	Acceptance Criteria	Example
Order of the peaks in the spectral profile (intensity vs pixel) from left to right	4-dye: blue-green-yellow-red	Elus Green Vellow Red
Extraneous peaks in the raw data profile (intensity vs scan)	None Note: The E5 profile may include extraneous peaks outside the matrix peak region, which can be ignored.	E5:
Peak morphology in the spectral profile (intensity vs pixel)	<ul> <li>No gross overlaps, dips, or other irregularities</li> <li>Peaks separate and distinct</li> <li>Peak apexes are separate and distinct (the tails will overlap)</li> </ul>	

- 3. (Optional) Review the CRL accuracy to determine discrepancies from the reference sequence:
  - General sequencing: 40 to 539 bp
  - MicroSeq<sup>®</sup> ID: 20 to 619 bp

If you observe large discrepancies (for example, 5 to 10 contiguous miscalled bases in the middle of a sequence), review the data. If you see a raw data peak larger than the adjacent peaks with baseline pull-up in all 4-dye color channels, it may indicate the presence of a bubble. Check the pump, run the Remove Bubbles wizard (see "Remove bubbles from the polymer pump" on page 251), then repeat the run as needed.

- 4. If the data for all capillaries meet the criteria above, click Accept Results.
- 5. If the data for the required number of capillaries do not meet the criteria above (7 capillaries for 8-capillary instruments, 21 capillaries for 24-capillary instruments):
  - a. (Optional) If you want to generate a report for the failed calibration, click **View Summary Report** or **View Detail Report** before you click Reject Results. To save the report electronically, select **CutePDF** as the printer.

b. Click **Reject Results**. For troubleshooting information, see "Sequencing install standard troubleshooting" on page 302.

**IMPORTANT!** If you reject results, the spectral calibration is not saved.



### Example sequencing install standard results

### View previously run sequencing install standards

Select **History View**, then select an install standard to view the associated calibration information.

### View and print a sequencing install standard report

**IMPORTANT!** Ensure that all dyes are selected before viewing the report. The report may contain incomplete data if all dyes are not selected.

Note the following:

• Install standard reports include the most recent install date if a capillary array was removed, then re-installed on the instrument. Spatial and spectral calibration reports include the date on which a capillary array is installed on the instrument for the first time.

- The sorting in the Install Standard screen is not applied to the report.
- You can generate a report for a failed installation standard run before you click Reject Results.
- 1. Click 📄 View Summary Report or 📄 View Detail Report.
- 2. In the Report screen, click toolbar options to manipulate the report as needed. Place the mouse pointer over an item for a description of the item.

Page 1 of 1	6 6 9 9 🗅 📄 🔊	:
	Modify settings for this report	
	🐸 Modify report settings 🛛 🛛	
:08-Feb-2	i one socialitys	
	Select the font to be used in reports.	
	10 🗸	
	OK Cancel	

- 3. To print the report, click 🌺 **Print**.
- 4. To save the report electronically (.pdf), print the report and select **CutePDF Writer** as the printer.
- 5. Close the report.



### Save historical performance check reports (.pdf) for record keeping

**IMPORTANT!** After performing a performance check, save the performance check report electronically for record keeping. The software does not save historical calibration results. Only the most recent spectral calibration for each dye set is maintained in the software.

- 1. Click 📄 View Summary Report or 📄 View Detail Report.
- 2. Click 🏊 Print.
- 3. In the Printer dialog box, select CutePDF Writer as the printer.
- 4. Specify a name and location for the report.

# Run the fragment analysis or HID Install standard performance check

### When to perform

When your instrument is installed, the service engineer runs a fragment analysis or HID install standard install performance check.

Applied Biosystems recommends that you run the fragment or HID install standard performance check monthly to verify that the instrument conforms to fragment analysis sizing precision, sizing range, and peak height specifications.

**IMPORTANT!** The performance check is application-specific. If you change polymer and capillary length, you must perform a new performance check.

Estimated run 30 minutes time

### Prepare for the fragment or HID install standard performance check

Prepare the instrument	<ol> <li>If you have not already done so, perform a spatial calibration (see "Spatial calibration" on page 99).</li> </ol>
	2. In the Dashboard, check consumable status (page 29). Ensure that:
	Consumables are not expired
	Adequate injections remain for consumables
	3. Ensure that the fluid levels are at the fill lines ("Check buffer fill levels" on page 31).
	4. Set the oven temperature to 60 °C, then click Start Pre-heat.
	Pre-heat the oven and detection cell while you prepare for a run (detection cell temperature is set by the software). Preheating helps mitigate subtle first-run migration rate effects. The pre-heat function automatically turns off after 2 hours.
	Applied Biosystems recommends that you pre-heat the oven for at least 30 minutes before you start a run if the instrument is cold.
	5. Check the pump assembly for bubbles and run the Remove Bubble wizard if needed (see page 251).
Prepare the installation standard plate	IMPORTANT! Do not use warped or damaged plates.

1. Prepare the installation standard as described in the product insert. See Table 29 on page 260 for standard part numbers.

Application	Installation Standard
Fragment analysis (G5 dye set, POP-7 <sup>™</sup> polymer, 50 cm capillary)	GeneScan <sup>™</sup> Installation Standard DS-33
HID (G5 dye set, POP-4 <sup>™</sup> polymer, 36 cm capillary)	AmpFtSTR <sup>®</sup> Identifiler <sup>®</sup> Allelic Ladder

2. Load the standards in injection position 1 in the plate:

**IMPORTANT!** You do not create a plate for the performance check. The software uses predetermined positions for the performance check run. You cannot specify standard location on the plate. If you do not place standards in the positions indicated, the calibration will fail.

8-capillary	A1 through H1
96-well plate	
24-capillary	A1 through H1, A2 through H2, and
96-well plate	A3 through H3
-	
	4
24-capillary	
384-well plate	and 5 in rows
•	A, C, E, G, I, K, N, O
Note: 384-well plates are not	
supported on	4
8-capillary	
instruments.	۵. ۱۳

- 3. Briefly centrifuge the plate containing the standards.
- 4. Verify that each sample is positioned correctly in the bottom of its well.
   IMPORTANT! If the reagents of any well contain bubbles or are not located at the bottom of the well briefly centrifuge the plate, remove the plate from the

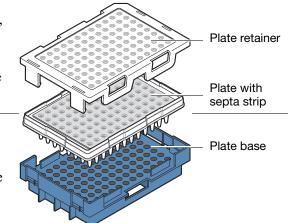
bottom of the well, briefly centrifuge the plate, remove the plate from the centrifuge, and verify that each sample is positioned correctly in the bottom of its well.

5. Store the plate on ice until you prepare the plate assembly and load the plate in the instrument.

Prepare the plate<br/>assemblyIMPORTANT! Prepare the plate assembly on a clean, level surface. Do not heat plates<br/>that are sealed with septa.

- 1. Align the holes in the septa strip with the wells of the plate, then firmly press downward onto the plate.
- 2. Align the holes in the septa strip with the wells of the plate, then firmly press downward onto the plate.
- 3. Place the sample plate into the plate base.

**IMPORTANT!** Make sure to use the correct plate base for standard plates versus 8-tube strips and fast plates. Using the wrong plate base may affect performance.



- 4. Snap the plate retainer (cover) onto the plate, septa, and plate base.
- 5. Verify that the holes of the plate retainer and the septa strip are aligned. If not aligned, re-assemble and then assemble the plate assembly.

**IMPORTANT!** The array tips will be damaged if the plate retainer and septa strip holes do not align correctly.

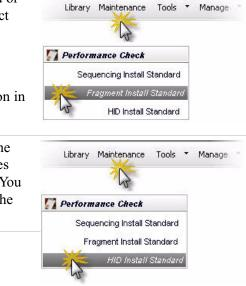
- Load the plate in the instrument
- 1. Place the plate in the autosampler with the labels facing you (or the instrument door) and the notched corner of the plate in the notched corner of the autosampler.
- AB 3500 96-Woll Fast 96-Woll Fast 96-Woll Fast 98-Woll Fast 98-Woll Fast 96-Woll Fast 96-Woll Fast 96-Woll Fast 86-00
- 2. Close the instrument door to re-initialize the instrument.

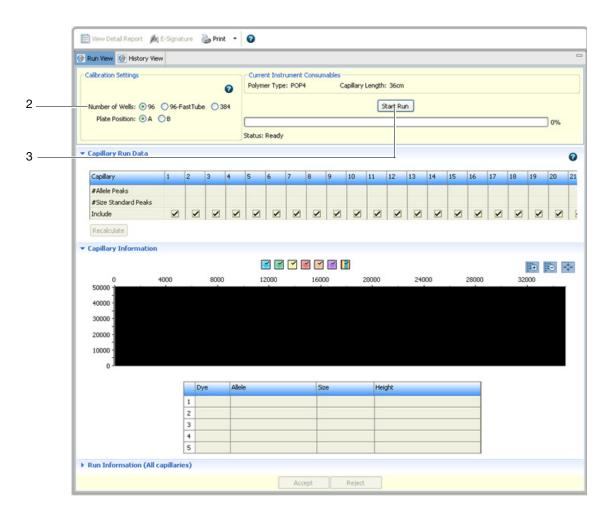
### Run the fragment analysis or HID install standard performance check

- Access the Fragment Install Standard or the HID install standard screen: Select Maintenance, then select Fragment Install Standard or HID Install Standard in the navigation pane.
- 2. Select the plate type and plate position in the instrument.

**Note:** You do not create a plate for the performance check. The software uses predetermined positions for the run. You cannot specify standard location on the plate.





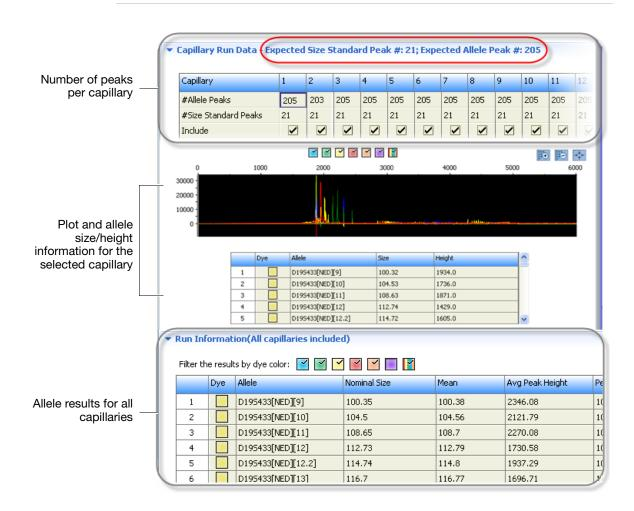


### What you see during a run

The system performs one run and indicates the number of observed allele and size standard peaks.

The Capillary Run Data display updates after the run is complete. The number of observed size standard and allele peaks is shown. Results for each allele are shown at the bottom of the screen in the Run Information table.

**Note:** The example shown below is for the HID install standard.



## How the software determines passing and failing capillaries for the fragment/HID performance check

The software evaluates peaks in the data for each capillary. To be identified as a possible allele, peaks must be within the following ranges (nominal allele size, or reference bin size, is hard-coded):

- All markers except THO1:  $\pm 0.7$  bp of nominal size for the allele
- THO1:  $\pm 0.5$  bp of nominal size for the allele

For all peaks that are within the nominal size range, the software calculates the Average Peak Height and the Sizing Precision. Peaks that meet the thresholds below pass.

Result	Description	Threshold		
Avg Peak Height	Average of peak heights for observed allele peaks of the included capillaries.	<ul> <li>Fragment: &gt; 175 RFU</li> <li>HID: &gt; 400 RFU</li> </ul>		
Sizing Precision	Standard deviation of the observed allele fragment sizes.	≤0.15 for expected alleles		
Pass/Fail	Alleles with a sizing precision and average peak height that do not meet thresholds fail <b>Note:</b> Review the data for failed alleles as described below.			

Result	Description				
For information only					
Nominal Size	Expected allele fragment peak size (bp).				
Mean	Average fragment size for the observed allele peaks.				
Peak Height % > Min	Percentage of observed allele peaks with a peak height above the minimum threshold.				
Sizing Accuracy	Difference between the allele size and the mean allele size.				

### Evaluate fragment/HID install standard data

1. Examine the number of size standard and allele peaks found for each capillary.

**Note:** The number of expected peaks shown below is for the HID install standard.

Expected –	- Capillary Run Data (E	xpected	Size S	itanda	rd Pea	<b>k #: 2</b> :	l; Expe	cted A	Allele P	'eak #	: 205	)	
	Capillary	1	2	3	4	5	6	7	8	9	10	11	12
Observed	#Allele Peaks	205	203	205	205	205	205	205	205	205	205	205	205
	#Size Standard Peaks	21	21	21	21	21	21	21	21	21	21	21	21
	Include	✓										-	V

2. If the expected number of alleles and size standard peaks are found, click Accept Results.

If the expected number of alleles and size standard peaks are not found, troubleshoot as described below.

Troubleshoot 1. Click a capillary with fewer than the Capillary expected number of #Allele Peaks 203 204 peaks to display #Size Standard Peaks 21 detailed information Include for each allele in the table below the plot. 2. Double-click the Size column to sort results and identify

> not found. A "0" Size value indicates that an allele falls outside the expected size window (Nominal Size  $\pm 0.7$  bp or  $\pm 0.5$  for THO1).

the alleles that were 5

Dye Allele Height D18551[NED][20] 0.0 1 0.0 2 D351358[VIC][17] 0.0 0.0 3 D195433[NED][9] 100.31 858.0 4 D195433[NED][10] 104.49 867.0 AMEL[PET][X] 105.01 931.0

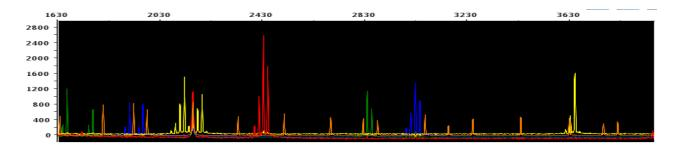
- 3. Troubleshoot failing data:
  - a. Analyze the install standard data files in your secondary analysis software (GeneMapper<sup>®</sup> Software v4.1 or later; GeneMapper<sup>®</sup> *ID-X* Software Software v1.2 or later using Identifiler<sup>®</sup> kit panels and bins).
  - b. Evaluate the failed data and examine the alleles not found by the 3500 Series Data Collection Software.
  - c. If the alleles are properly called in the secondary analysis software, you can:
    - Deselect the Include checkmark for a capillary.
    - Click Recalculate.
    - Accept the install standard results.

Note: The GeneMapper<sup>®</sup> *ID-X* Software may identify alleles not identified by the 3500 Series Data Collection Software because of the bin-offsetting feature (which uses the observed alleles in the allelic ladder samples to adjust the reference bin locations for samples).

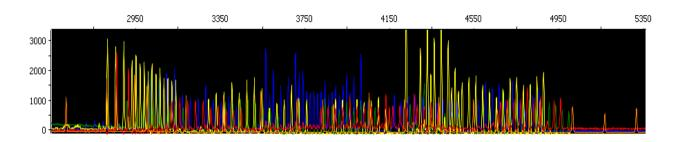
If the alleles are not properly called:

- (Optional) Click View Detail Report to save a record of the failed run. To save the report electronically (.pdf), print the report and select CutePDF Writer as the printer. For more information, see "Save historical performance check reports (.pdf) for record keeping" on page 138.
- Click Reject Results.
- Rerun the install standard to determine if the problem may be caused by sample preparation, a poor injection, a capillary issue, or a system problem (which may require instrument service). For more information, see "Fragment/HID install standard troubleshooting" on page 303.

### Example fragment install standard results



### Example HID install standard results



### View previously run install standards

Select **History View**, then select an install standard to view the associated calibration information.

### View and print a fragment or HID install standard detail report

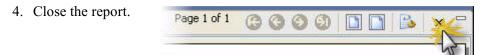
**IMPORTANT!** Ensure that all dyes are selected before viewing the report. The report will contain incomplete data if all dyes are not selected.

Note the following:

- Install standard reports include the most recent install date if a capillary array was removed, then re-installed on the instrument. Spatial and spectral calibration reports include the date on which a capillary array is installed on the instrument for the first time.
- The sorting in the Install Standard screen is not applied to the report.
- To generate a report for a failed installation standard run, you must do so before you click Reject Results.
- 1. Click 📄 View Detail Report.
- 2. In the Report screen, click toolbar options to manipulate the report as needed. Place the mouse pointer over an item for a description of the item.



3. To print the report, click 🌺 **Print**.



### Save historical performance check reports (.pdf) for record keeping

**IMPORTANT!** After performing a performance check, save the performance check report electronically for record keeping. The software does not save historical calibration results. Only the most recent spectral calibration for each dye set is maintained in the software.

- 1. Click 🔛 View Detail Report.
- 2. Click 🏊 Print.
- 3. In the Printer dialog box, select CutePDF Writer as the printer.
- 4. Specify a name and location for the report.

### **Overview of libraries**

The Library workflow contains the screens where you manage assays, protocols, and other items that you use to acquire and process data.

The Library workflow contains:

- Items that you select when you set up a run:
  - Plates
  - Assays
  - Optional filename conventions
  - Optional results groups
- Items that you select when you create an assay:
  - Instrument protocols
  - Primary analysis protocols Basecalling (sequencing), sizecalling (fragment analysis), QC (HID analysis)
  - Optional secondary analysis protocols Sequencing analysis, fragment analysis, and HID analysis
- Items you select when you create instrument sizecalling and QC protocols:
  - Dye sets
  - Size standards

#### Factory-provided, template, and locked items

The 3500 Series Data Collection Software libraries include factory-provided items that are optimized for different applications (for example, instrument protocols with specific run modules and primary analysis protocols with specific settings). You can use the factory-provided items directly. If the factory-provided items do not suit your needs, you can modify the factory-provided items, or create new items.

Entries in the library may be flagged with the following symbols:

- Factory-provided. Cannot be edited or deleted.
- 📝 Template.
- Locked. If your system includes the SAE module, can be unlocked and modified only by the user who created it, the administrator, or another user with unlock permissions. For information, see Chapter 7, "Use Security, Audit, and E-Sig Functions (SAE Module)" on page 197.

### General library procedures

### **Access libraries**

Select **Library** in the menu bar to access the Library workflow.



The Library workflow contains the screens where you manage assays, protocols, and other items that you use to acquire and process data.

The Library workflow contains:

- Items that you select when you set up for a run: plates, assays, filename conventions, and results groups
- Items that you select when you create an assay:
  - Instrument protocols
  - Primary analysis protocols Basecalling (sequencing), sizecalling (fragment analysis), QC (HID analysis)



- Optional secondary analysis protocols –
   Sequencing analysis, fragment analysis, and HID analysis
- Items you select when you create instrument, sizecalling, and QC protocols:
  - Dye sets
  - Size standards

You can click **Main Workflow**, or select **Dashboard** or any other menu item at any time to advance from the Library workflow.

### Create a new entry from a factory-provided, template, or locked entry

**IMPORTANT!** Auditing of an item depends on whether it is created directly from the library or from within another item (for example, you can create an assay directly from the library, or within a plate in the Assign Plate Contents screen). For more information on auditing, see "Review the object audit history" on page 210.

- 1. Select the factory-provided entry in the library.
- 2. Click W Duplicate. The software creates a "Copy of" the item you duplicated.
- 3. Select the "Copy of" item, then click 🥖 Edit.
- 4. Enter a name for the item.
- 5. Modify parameters as needed (see the appropriate section for information).

6. Click Save.

#### Delete a library entry

**IMPORTANT!** Auditing of an item depends on whether it is deleted directly from the library or from within another item (for example, you can delete an assay directly from the library, or within a plate in the Assign Plate Contents screen). For more information on auditing, see "Review the object audit history" on page 210.

**Note:** You cannot delete states factory-provided items.

Select an item, then click 🕮 Delete.

Deleting a library entry does not affect existing items that contain the entry. (When you select an item to include in a higher-level item, a *copy* of that item is included in the higher-level item. For example, when you select an instrument protocol to include in an assay, a copy of the instrument protocol is included in the assay. If you delete the instrument protocol, the copy of the instrument protocol in the assay remains intact.)

For information on how deleted items are tracked in auditing, see "Audit action" on page 210.

### Edit a library entry

**IMPORTANT!** Auditing of an item depends on whether it is edited directly from the library or from within another item (for example, you can edit an assay directly from the library, or within a plate in the Assign Plate Contents screen). For more information on auditing, see "Review the object audit history" on page 210.

- 1. Select an item, then click 🌽 Edit.
- 2. Modify parameters as needed.
- 3. Click Save.

### Import and export a library entry

You can import and export .xml files for use with other 3500 or 3500xL analyzer instruments:

- **Import** Click **Import**, then select the .xml file to import. If any items in the import file exist in the library, the software displays a message and gives you the option to replace or skip the item.
- **Export** Select one or more entries, then click **Export**, then specify a location for the export file.

To select multiple entries, Shift-click to select contiguous entries, Ctrl-click to select non-contiguous entries.

### View audit and e-signature histories for library entries

**Note:** An administrator can also view audit and e-signature histories in the SAE module. For information, see Chapter 7, "Use Security, Audit, and E-Sig Functions (SAE Module)" on page 197.

To view the audit or e-signature history for a library entry:

- 1. Select the item in the library.
- 2. Click View Audit History or View E-Signature History (active only if the selected item is enabled for e-sig).

**Note:** Factory-provided items do not list creation date in the audit history. If you duplicate a factory-provided item, the new item contains an audit history that starts with the duplication date listed as the creation date.

3. For more information, see "Display audit histories" on page 209.

#### Sort, filter, and search library entries

**Sort** Double-click column headers to sort. Multi-column sorting is supported:

- Double-click a column header to sort the column.
- Alt+Shift-click another column header to sort another column.
- Alt+Shift-click a third column header to sort a third column.

Numbers in the column headers reflect sort order.





You can select an application type from the Filter list to display only plates for the selected application.

Search

Search:		All	¥
		All	~
		Plate Name	
		Type Description	
Desc	ription	Description	_
	·	Barcode	×

In each library, you can select a category to search, then enter the text to search for. The list of categories corresponds to the column headers in each library.

Click **Go** to search. Click **Clear** to remove the search criteria.

### Customize a library table

Click the Table Settings button, then specify the columns to show or hide.

Click:

Save to

• Apply – To use the settings for this session only.

Column A	Column B	Calumn C	Column L
Table Preference	es		
Select your table display	preferences below.		
Available Columns to	Display		
Column A	<u>^</u>		
Column B			
Column C		Add Select	ed >>
Column D		Add All	The second se

Preferences – To

save for future use by all users. If your system includes the SAE module, preferences are saved for the logged-in user.

• Restore Defaults – To restore factory default settings.

### **Plates library**

The Plates library contains all plates that have been saved in the software (plates that have been run and plates that have not yet been run).

### **Plate overview**

**Plate definition** A plate associates sample attributes (sample information and analysis information) with a well position. A plate defines how samples are analyzed during capillary electrophoresis and how sample files are named and stored after analysis.

When you create a plate, you specify:

- Plate type (sequencing, fragment, mixed, or HID)
- Number of wells, capillary length, and polymer type

When you set up a plate for a run, you add assays, optional file name conventions, and optional results groups to wells in the plate. If you add these items from the library, a *copy* of the items is added to the plate, and can be modified independently from the original items stored in the library. For information on how changes are tracked if auditing is enabled, see "Audit action" on page 210.

**Plate templates** The Plates library includes templates that are optimized for different applications (for example, plates defined with the appropriate polymer and capillary length) that you can use to create new plates.

### Create a new plate

- 1. Access the Plates library.
- 2. Click 🚟 Create.

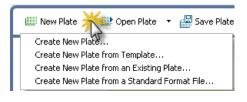
The software switches to the Main workflow and displays the Define Plate Properties screen (Figure 11 on page 145).

**Note:** You can also access the Define Plate Properties screen from the Dashboard and the Assign Plate Properties screen.



3. To create a new plate, specify settings (Table 6 on page 145).

To create a new plate based on an existing plate, click **New Plate**, then select an option. Select a plate, click **Open**, then specify settings.



For information on other Create New Plate options, see:

- "Create a plate from a template" on page 43.
- "Create a plate for importing" on page 73 (for the Create New Plate from a Standard Format File option)
- 4. Select a Save option.



📖 New Plate 🔹 🔜 Open Plate 🔹 🔛 Save Plate 🔹 📑 Close Plate	🛃 Start Run
Plate Details	0
* Name: * Number of Wells: Plate Name is a required field. Provide a unique value. * Plate Type: HID * Capillary Length: 36 * Polymer: POP4 * Polymer: POP4	Owner:Barcode:
▼ Secondary Analysis	Perform Auto-Analysis (
* Software Type:	
* Username:	
* Password:	
* Auto Analysis is performed: Only when the results group is complete (	When every injection completes

Figure 11 Define Plate Properties

#### Table 6Define Plate Properties

Setting	Description			
Plate Details				
Name	Plate name. Names must be unique.			
Number of Wells	• 96 well – For standard 96-well plates			
	<ul> <li>96–Supports 96-well standard reaction plate. 8-strip standard tubes are also supported with appropriate retainers.</li> </ul>			
	• 96 Fast tube – For Fast 96-well plates and 8-strip tubes			
	<ul> <li>96-Fast Tube– Supports 96-well Fast reaction plate. 8-strip fast tubes are also supported with appropriate retainers.</li> </ul>			
	• 384 well – 384-well plates			
Plate Type	Sequencing			
	Fragment analysis			
	Mixed (Seq + Frag)			
	• HID			
Capillary Length and Polymer	Capillary length and polymer type with which the plate will be used			
Owner, Barcode, Description (optional)	Optional text entries			
	You can use these entries to search for plates in the Plates library and in run logs (Tools ► View Logs).			

#### Table 6 Define Plate Properties (continued)

Setting	Description

#### Secondary Analysis

**Note:** The secondary analysis protocol settings you specify in an assay must match the auto-analysis settings for the plate. For more information, see:

- "Sequencing analysis protocols library (secondary analysis)" on page 189
- "MicroSeq<sup>®</sup> ID protocols library (secondary analysis)" on page 191
- "Fragment analysis protocols library (secondary analysis)" on page 193
- "HID analysis protocols library (secondary analysis)" on page 195

Perform Auto-Analysis	Enables the plate for use with auto-analysis with a supported secondary software
Software Type	Supported secondary software
Software Location	Computer on which the supported secondary software is installed
Username	User name and password required by the secondary analysis software
Password	
Auto-Analysis is performed	Determines when data is sent to the secondary analysis software:
(fragment/HID only)	Only when the results group is complete
	When every injection completes

### **Assays library**

### Assay overview

An assay contains the instrument protocol (dye set and run configuration) and primary analysis protocol needed to collect data and basecall or sizecall a sample. Assays, File Name Conventions, and Results Groups may already be listed in the plate template when you create a plate from a template.

Note: If no assay is listed, add at least one assay.

An assay contains:

- One or more instrument protocols appropriate for the sample type/dye set for which the assay will be used
- A primary analysis protocol that depends on your application:
  - Sequencing Basecalling protocol
  - **Fragment** Sizecalling protocol
  - HID QC protocol
- (Optional) A secondary analysis protocol that depends on your application:
  - Sequencing SeqScape<sup>®</sup> Software v2.7 or later) or MicroSeq<sup>®</sup> ID Analysis Software v2.2 (or later)
  - **Fragment analysis** GeneMapper<sup>®</sup> Software v4.1 (or later)
  - **HID** GeneMapper<sup>®</sup> *ID-X* Software ID-X Software v1.2 (or later)

Assays are required by all application types. You must assign an assay to all named sample wells on a plate before you can link a plate and run it.

When you create an assay, you add one or more instrument protocols and a primary analysis protocol. If you add these items from the library, a *copy* of the items is added to the assay, and can be modified independently from the original items stored in the library. For information on how changes are tracked if auditing is enabled, see "Audit action" on page 210.

### Create a new assay

If factory-provided assays do not suit your needs, you can create new assays:

- 1. Access the Assays library. 블 3500 Data Collection Software 2. Click **The Create**. Dashboard Edit 🔻 Library Maintenan 🚽 🕁 eate Note: You can also create an Library Resources assay from the Assign Plate Contents screen. Filter: All 🕻 Manage Plates Assay Name 1 AS YF\_POP4 File Name Conventions 100 New Plate + 100 Open Plate + 200 Save Plate + 100 Open Plate To state inere EE Table Vie Setup 212 Define Plate Properties Assign Plate Content Her Assays Actions \* New 55+Normaliz... h Edit Duplicate Rename Delete Add From Library Save To Library Import From File Select Related Wells
- 3. In the Create New Assays dialog box, select an application type: Sequencing, Fragment, or HID. The screen changes depending on the application type you select (Figure 12 on page 149 shows the sequencing screen).
- 4. Specify settings (see Table 7 on page 149).
- 5. Save the assay:
  - If you are creating the assay from the Library, click Save.
  - If you are creating the assay from the Assign Plate Contents screen, click **Apply to Plate** or **Save to Library**.

🔚 Create New Assay					
Setup an Assay ⑧ Basecalling Protocol cannot be empty.					
* Assay Name: New Assay Application Type: Sequencing V		Locked	Color:	Black	<ul> <li></li> </ul>
Do you wish to assign multiple instrument protocols	to this assay? 💿 No	🔘 Yes			
* Instrument Protocol:			*	Edit	Create New
* Basecalling Protocol:			*	Edit	Create New
SeqScape/MicroSeqID Protocol:			¥	Edit	Create New
Close					Save

Figure 12 Create New Assay – sequencing – the highlighted area changes based on the Application Type

#### Table 7 Assay settings

Setting	Description						
Assay Name	Name of the assay. Names must be unique.						
Locked	When enabled, allows the entry to be unlock it, the administrator, or another user with unl includes the SAE module (described in Chap Functions (SAE Module)" on page 197.	ock	ре	rmissions. Us	eful when	your sys	
Color	Color code for the assay when it is		1	2	3	4	5
	displayed in the Assign Plate Contents screen (if Assay Color is selected for Show In Wells).	A	H	H	H	Н	H
		в	H	н	н	н	B
		-	н		Assa	ys	
		c					Actions
				🖲 🗌 🗛 HID:	36_POP4_G5+	No 📝	23
		15	ан	🖲 🗌 🛤 HID:	36_POP4_G5_	ls 📝	13
				and the second s	36_POP4_G5_	and the second second	
				🕘 🗌 🙈 HID:	36_POP4_G5+	No 🥖	13
Application Type	Sequencing						
	Fragment analysis						
	• HID						

### Table 7 Assay settings (continued)

Setting		Description		
Do you wish to assign multiple instrument protocols to this assay?	When you select Yes, allows you to select or create additional instrument protocols for the assay. The software creates one injection for each instrument protocol specified in an assay.	Protocols         Do you wish to assign multiple instrument protocols to this assay?         Instrument Protocols         Available Library         Instrument Protocols:         O Instrument Protocols:         O Instrument Protocol(s) Assigned to this Assay         Edit       Remove         Move Up       Move Down         NOTE: Order the list of protocols in the order you want them injected		
Instrument Protocol	Instrument protocol fo	r data collection.		
	For information, see "I	nstrument protocol library" on page 165.		
Sequencing				
Basecalling Protocol	Protocol for primary analysis (basecalling and trimming) and quality determination.			
	For information, see "E page 174.	Basecalling protocols library (primary analysis – sequencing)" on		
SeqScape software	Optional protocol for s	econdary analysis (auto-analysis).		
/MicroSeq <sup>®</sup> software D Protocols	For information, see:			
		sis protocols library (secondary analysis)" on page 189.		
	"MicroSeq <sup>®</sup> ID prot	ocols library (secondary analysis)" on page 191.		
Fragment				
<ul> <li>Sizecalling Protocol</li> </ul>	Protocol for primary ar	nalysis (peak detection and sizing) and quality determination.		
	For information, see "Spage 179.	Sizecalling protocols library (primary analysis – fragment)" on		
GeneMapper <sup>®</sup>	Optional protocol for s	econdary analysis (auto-analysis).		
software Protocol	For information, see "F page 193.	Fragment analysis protocols library (secondary analysis)" on		
HID				
QC Protocol	Protocol for primary ar	nalysis (peak detection and sizing) and quality determination.		
	For information, see "	QC protocols library (primary analysis – HID)" on page 184.		
GeneMapper <sup>®</sup> ID-X	Optional protocol for s	econdary analysis (auto-analysis).		
Protocol For information.		HD analysis protocols library (secondary analysis)" on page 195.		

### File name conventions library

### File name convention overview

A File Name Convention (FNC) specifies the naming convention for sample data files. It is an optional component in a plate.

If you do not specify a file name convention, data files are named in this format:

<sample name>\_<well>

The file extension is determined by the application you run:

- Sequencing .ab1 (you can also set Preferences to export additional file formats. See "Set sequencing preferences" on page 36.)
- Fragment analysis .fsa
- **HID** .hid

**Note:** The file location specified in a file name convention is used only if a results group is not specified for a well.

When you set up a plate for a run, you can optionally add file name conventions to the plate. If you add this item from the library, a *copy* of the item is added to the plate, and can be modified independently from the original item stored in the library. For information on how changes are tracked if auditing is enabled, see "Audit action" on page 210.

### Create a new file name convention

If factory-provided file name conventions do not suit your needs, you can create new file name conventions:

- 1. Access the File Name Conventions library.
- 2. Click 式 Create.

**Note:** You can also create a file name convention from the Assign Plate Contents screen.



Districted	10 New Place + 12 Open Place + 20 Law Place + 12 Open Place	
	Important         Important <t< th=""><th>ions Actions 🔹</th></t<>	ions Actions 🔹
- Contraction of the second se	eneral File Name	New Edit Duplicate Rename Delete Add From Library Save To Library Import From File

3. In the Create New File Name Conventions dialog box (Figure 13 on page 153), select attributes and delimeters (see Table 8 on page 153).

As you select attributes, the software displays a preview of the file name.

Available Attributes	- Selected Attributes	
Amplicon Name Analysis Protocol Name Assay Name Capillary Number Custom Text1 Custom Text2	Add >>> Sample Name	
Custom Text3 Date of Run	Move Up	
Injection Number Instrument Name Instrument Protocol	Move Down	
Owner Name - Select File Name	Attributes	
Preview of File	Name: <pre>Sample Name&gt;<capillary number=""></capillary></pre>	
-Available Attr	ibutes	Selected Attributes

- 4. To add delimiters between items in the Selected Attributes list:
  - a. Ctrl-click or Shift-click to select two or more attributes.
  - b. Select a delimiter.
  - c. Select the Add between attributes check box.
  - d. Click Add.
- 5. Save the file name convention:
  - If you are creating the file name convention from the Library, click Save.
  - If you are creating the assay from the Assign Plate Contents screen, click **Apply to Plate** or **Save to Library**.

Create New File Name Convention
Setup a File Name Convention       Image: Setup a File Name Convention         Name is a required field. Provide a unique value.       Image: Setup a File Name Convention
* Name: Locked Color: Black
Select File Name Attributes
Preview of File Name: <sample name=""></sample>
Available Attributes   Amplicon Name   Analysis Protocol Name   Assay Name   Capillary Number   Custom Text1   Custom Text2   Custom Text3   Date of Run   Injection Number   Instrument Protocol   Owner Name   V Delimiters   Select a delimiter   Plus ( + )   Add >>
Add a custom value to available attributes (optional) Custom Text 1: Custom Text 2: Custom Text 3:
Select File Location C:\Applied Biosystems\3500\Data
Custom File Location Browse
Close
Close

Figure 13 Create New File Name Convention

Table 8	File name	conventions	settings
---------	-----------	-------------	----------

Setting	Description
Name	Name of the file name convention. Names must be unique.
Locked	When enabled, allows the entry to be unlocked and modified only by the user who created it, the administrator, or another user with unlock permissions. Useful when your system includes the SAE module (described in Chapter 7, "Use Security, Audit, and E-Sig Functions (SAE Module)" on page 197.

Setting	Description
Color	Color code for the file name convention when it is displayed in the Assign Plate Contents screen (if File Name Convention Color is selected for Show In Wells).
Preview of name Available attributes	Interactively displays the attributes you select.         • Amplicon Name (from Customize       • Polymer Type
	<ul> <li>Sample Info in sequencing assays)</li> <li>Analysis Protocol Name – (primary analysis protocol)</li> <li>Assay Name</li> <li>Capillary Number</li> <li>Custom Text fields (up to 3)</li> <li>Date of Run</li> <li>Injection Number</li> <li>Instrument Name</li> <li>Instrument Protocol</li> <li>Owner Name (plate owner)</li> <li>Plate Name</li> <li>IMPORTANT! The maximum allowed length of a file name, including the path, is 240 characters. The software warns you if your selections will possibly exceed the maximum, but allows you to save the file name convention. However, you will see a pre-check validation error when you start a run if the file name will exceed 240 characters.</li> </ul>
Delimiters	Symbols you can include in the file name: Dash (-), Dot (.), Underscore (_), Plus (+), Dollar (\$).
Custom text	Text to display for the custom text attribute fields.
File location	<ul> <li>The file location in the file name convention is used only if no results group is specified for a well.f</li> <li>The Results Group file location overrides the File Name Convention file location.</li> </ul>

### Table 8 File name conventions settings (continued)

### **Result group library**

### **Results group overview**

A Results Group is used to name, sort, and customize the folders in which sample data files are stored. It is an optional component in a plate.

**Note:** The file location specified in a results group overrides the file location in the file name convention specified for a well.

When you set up a plate for a run, you can optionally add results groups to wells in the plate. If you add this item from the library, a *copy* of the item is added to the plate, and can be modified independently from the original item stored in the library. For information on how changes are tracked if auditing is enabled, see "Audit action" on page 210.

#### Allelic ladder location (HID analysis)

To accurately genotype samples, the GeneMapper<sup>®</sup> ID-X Software requires at least one allelic ladder sample per run folder. (Multiple allelic ladder samples in a single run folder can also be used for analysis.)

Applied Biosystems recommends that you run one allelic ladder for 24 a set of samples:

- 8-capillary instruments One allelic ladder per 3 injections
- 24-capillary instruments One allelic ladder per 1 injection

**Note:** Run HID validation studies to determine the required number of allelic ladders for your application.

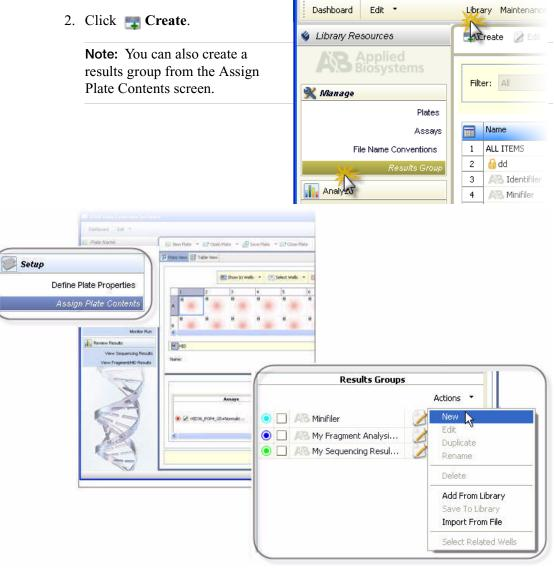
See "Results group example 2: store one allelic ladder per run folder (8-capillary instruments)" on page 161 for a results group example that places three injections in each run folder for 8-capillary instruments.

### Create a new results group

If the factory-provided results groups do not suit your needs, you can create new results groups:

블 3500 Data Collection Software

1. Access the Results Groups library.



3. In the Create Results group dialog box (Figure 14 on page 158), select attributes and delimeters (see Table 9 on page 158).

As you select attributes, the software displays a preview of the results group name.

Available Attributes PA Protocol Name Plate Name Prefix Delimiters Celect Results Group Attributes	Add >> Celected Attributes Results Group Name
Preview of Results Group Name	e: <results group="" name=""><pa name="" protocol=""><plate name=""></plate></pa></results>
Available Attributes Prefix	Selected Attributes Results Group Name

- 4. To add delimiters between items in the Selected Attributes list:
  - a. Ctrl-click or Shift-click to select two or more attributes.
  - b. Select a delimiter.
  - c. Select the Add between attributes check box.
  - d. Click Add.
- 5. Save the results group:
  - If you are creating the results group from the Library, click Save.
  - If you are creating the results group from the Assign Plate Contents screen, click **Apply to Plate** or **Save to Library**.

The Results Group file location overrides the File Name Convention file location.

🔄 Create New Results Group
Setup a Results Group       Image: Setup a Results Group         Image: Name is a required field. Provide a unique value.       Image: Setup a Results Group
0
* Name: Locked Color: Black
Select Results Group Attributes
Preview of Results Group Name: <results group="" name=""></results>
Available Attributes       Add >>       Selected Attributes         Assay Name       Add >>       Results Group Name         Ipection Number       <
Select a delimiter Dash ( - ) Move Down Add between attributes Add >>
Enter a custom value as either the Prefix or Suffix (optional) Prefix: Suffix:
Select Reinjection Folder Option
⊙ Store reinjection sample files in a separate Reinjection folder (same level as Injection folders)
◯ Store reinjection sample files with original sample files (same level)
C Select Folder Option
Default file location C:\Applied Biosystems\3500\Data\ <ir folder="">\<results folder="" group="" name="">\<inj folder="">\</inj></results></ir>
O Custom file location
🗹 Include an Instrument Run Name folder
☑ Include a Result Group Name folder
Include an Injection folder
Close Save

Figure 14 Create New Results Group

#### Table 9 Results group settings

Setting	Description
Name	Name of the results group. Names must be unique.
	The Results Group Name is a required attribute, you cannot remove this attribute from the Selected Attribute list.
Locked	When enabled, allows the entry to be unlocked and modified only by the user who created it, the administrator, or another user with unlock permissions. Useful when your system includes the SAE module (described in Chapter 7, "Use Security, Audit, and E-Sig Functions (SAE Module)" on page 197.

Setting	Description					
Color	Color code for the results group when it is displayed in the Assign Plate Contents screen (if Results Group Color is selected for Show In Wells).	1         2         3           A         B         B           C         Results Groups			4	
		D	<ul> <li>AS Minifiler</li> <li>AS My Fragm</li> <li>AS My Seque</li> </ul>	100	Action	
Preview of name	Interactively displays the attributes you select.					
Available attributes	<ul> <li>Results Group Name (required)</li> <li>Assay Name</li> <li>Injection Number</li> <li>IP Name (instrument protocol)</li> <li>Logged-in User Name (available only when security is enabled in the SAE module)</li> </ul>	•   •   • {	PA Protocol Name ( Plate Name Prefix Start Instrument Ru Suffix		-	
Delimiters	Symbols you can include in the results group name: Dash (-), Dot (.), Underscore (_), Plus (+), Dollar (\$).					
Prefix/suffix text	Text to display for the prefix or suffix text attribute fields.					
Select re-injection folder option	<ul> <li>Store reinjection sample files in a separate reinjection folder (same level as injection folders)</li> <li>Store reinjection sample files with original sample files (same level)</li> </ul>					
Select folder option	<ul> <li>Location:</li> <li>Default file location (specified in Preference)</li> <li>Custom location</li> <li>Sub-folder options:</li> <li>Include an instrument run name folder for Run screen)</li> <li>Include a results group name folder</li> </ul>			ned in the Loa	d Plates	
	<ul> <li>Include an injection folder</li> </ul>					

### Table 9 Results group settings (continued)

### Results group example 1: store files by plate name

Two default, factory-provided, results groups are provided that store sample data files by plate name:

- Figure 15 on page 160 shows the factory-provided PN\_Injfolder\_RG results group and the folders created when it is used. This results group creates a folder for each injection.
- Figure 16 on page 161 shows the factory-provided PN\_RG results group and the folders created when it is used. This results group does not create a folder for each injection. All samples for a plate are stored in the same folder. If you include two plates in a run, a separate folder is created for each plate.

-	
Edit Results Group PN_Injfolder_NG	s 🗀 C:\Applied Biosystems\3500\Data\Plate01-PN_Injfolder_RG\Inj1 2009-03-22-13-19-21-211
Setup a Results Group         * Name:       PN_Injfolder_RG         * Name:       PN_Injfolder_RG         * Name:       PN_Injfolder_RG         * New of Results Group Name       Selected Attributes         * Prever of Results Group Name       Plate Name         * Default file       Plate Name         * Default file       Selected Attributes         * Results Group Name       Plate Name         * Default file       Nor Use         * Store reinjection sample files with original sample files (same let         * Default file location       C:\Applied Biosystems\3500\Data\ <plate< td="">         * Include an Instrument Run Name folder       ************************************</plate<>	<ul> <li>Applied Biosystems</li> <li>3500</li> <li>Data</li> <li>Inji 2009-03-22-13-19-21-211</li> <li>Inji 2009-03-22-13-19-21-227</li> <li>Inji 2009-03-22-13-19-21-243</li> <li>Inji 2009-03-22-13-19-21-243</li> <li>Inji 2009-03-22-13-19-21-243</li> <li>Inji 2009-03-22-13-19-21-243</li> <li>test_D01.hid</li> <li>test_D01.hid</li> <li>test_D01.hid</li> <li>test_D03.hid</li> </ul>

Figure 15 PN\_Injfolder\_RG results group

Edit Results Group PN. Injfolder. RG		
Setup a Results Group	C:\Applied Biosystems\3500\Data\Plate02-PN_RG	
		× 🖻 AL 01_A01.hid
Name: PN_RG     Identify Selected Attributes     Plate Name     Plate Name     Plate Name     Plate Name     Dash ( - )     Results Group Name     Selected Attributes     Plate Name     Dash ( - )     Results Group Name     Selected Attributes     Plate Name     Dash ( - )     Results Group Name     Selected Attributes     Plate Name     Dash ( - )     Results Group Name     Selected Attributes     Plate Name     Dash ( - )     Results Group Name     Selected Attributes     Plate Name     Dash ( - )     Results Group Name     Selected Attributes     Plate Name     Dash ( - )     Results Group Name     Selected Attributes     Plate Name     Dash ( - )     Results Group Name     Dash ( - )     Results Group Name     Dash ( - )     Include an Instrument Run Name folder     Include an Instrument Run Name folder     Include an Injection folder	<ul> <li>Applied Biosystems</li> <li>3500</li> <li>Data</li> <li>Plate02-PN_RG</li> <li>Plate03-PN_RG</li> </ul>	<ul> <li>Interst_A02.hid</li> <li>Itest_A03.hid</li> <li>Itest_B01.hid</li> <li>Itest_B02.hid</li> <li>Itest_B03.hid</li> <li>Itest_B03.hid</li> <li>Itest_C01.hid</li> <li>Itest_C02.hid</li> <li>Itest_C03.hid</li> <li>Itest_C03.hid</li> <li>Itest_C03.hid</li> <li>Itest_C03.hid</li> <li>Itest_D03.hid</li> </ul>

Figure 16 PN\_RG results group

## Results group example 2: store one allelic ladder per run folder (8-capillary instruments)

Applied Biosystems recommends that you run one allelic ladder for each set of 24 samples (see "Allelic ladder location (HID analysis)" on page 155).

To store one allelic ladder per run folder on an 8-capillary instrument, create one results group for each set of three injections on the plate. Each results group specifies a results group name folder. Because you assign one results group to a set of three injections, all 24 sample data files, including the allelic ladder, are stored in the same results group folder.

The example below shows one results group; for a full 96-well plate, create three more with the same settings, but different names, for example, Injection 4 - 6, Injection 7 - 9, and Injection 10 -12.

🕒 Create	New Results Group					
* Name:	Injection 1 - 3	Locked				
-Select Re	Select Results Group Attributes					
Preview	Preview of Results Group Name: Injection 1 - 3					
Available Attributes       Add >>       Selected Attributes         IP Name       Add >>       Results Group Name         PA Protorol Name       <						
C Select Folder Option						
Operault file location C:\Applied Biosystems\3500\Data\ <ir folder="">\Injection 1 - 3\</ir>						
Custom file location						
✓ Include an Instrument Run Name folder						
☑ Include a Result Group Name folder						
Include an Injection folder						
Close		Save				

### Results group example 3: store re-injections in separate folders

Figure 17 on page 163 shows an example results group that specifies a sample file storage location of:

C:\Example\instrument run (IR) folder\result group name folder [results group name+start instrument run date/time stamp+logged in user name]\injection name or re-injection name folder.

The numbers in the figure relate the elements in the results group with the elements in the file hierarchy created by a run that uses this results group (Figure 20 on page 164).

* Name: RG Example	3
Selected Attributes Results Group Name Plus ( + ) Start Instrument Run Date/Time Stamp Plus ( + ) Logged in User Name	3
Select Reinjection Folder Option <ul> <li>Store reinjection sample files in a separa</li> <li>Store reinjection sample files with origin</li> </ul>	ate Reinjection folder (same level as Injection folders)
<ul> <li>Select Folder Option</li> <li>Default file location C:\Applied Biosyst</li> <li>Custom file location C:\Example</li> </ul>	ems\3500\Data\
<ul> <li>Include an Instrument Run Name folder</li> <li>Include a Result Group Name folder</li> <li>Include an Injection folder</li> </ul>	2 3 4

#### Figure 17 Results group example

Figure 18 on page 163 shows the injection list for a run that specifies duplicate and re-injections.

The numbers in the figure relate the elements in the injection list with the elements in the file hierarchy created by this run (Figure 20 on page 164).

onr	nection Statu Run Nam		nected 1 2009-02-05-14-59-56-7	703 5			Name: Adminis Itatus: Running	
	njection Lis jections crea		ails ' in Plate A - 0 in Plate B	(name used Instrument R Name folde	un			(logged-in user name)
F	Injection	Туре	Assay		Instrument Protocol	Plate	Analysis	Flags
ł	(duplicate and re-		IF+Norm_POP4_xl		HID36_POP4xl_G5	Plate 01	✓	
2	injections)		IF+Norm_POP4_xl		HID36_POP4×l_G5	Plate 01	✓	
3	<b>~</b>	₩	IF+Norm_POP4_xl		HID36_POP4×l_G5	Plate 01	✓	
4		₽.	IF+Norm_POP4_xl		HID36_POP4×l_G5	Plate 01	✓	
5	نے ا	₽.	IF+Norm_POP4_xl		HID36_POP4×l_G5	Plate 01	✓	
6	✓		IF+Norm_POP4_xl		HID36_POP4xl_G5	Plate 01	-	
7			IF+Norm_POP4_xl		HID36_POP4xl_G5	Plate 01		

#### Figure 18 Injection list example

Figure 19 on page 164 shows an example file name convention that specifies a sample name syntax of:

*sample name.(primary) analysis protocol name.unique time stamp integer* 

The numbers in the figure relate the elements in the file name convention with the files created by a run that uses file name convention (Figure 20 on page 164).

* Name: FNC Example	•		Locked
Select File Name Attribu	tes		
Preview of File Name:	<sample name="">.<ana< th=""><th>lysis Protocol Name</th><th>&gt;.<unique integer="" stamp="" time=""></unique></th></ana<></sample>	lysis Protocol Name	>. <unique integer="" stamp="" time=""></unique>
Available Attributes – Amplicon Name Assay Name Capillary Number Custom Text1 Custom Text2 Custom Text3		Add >> << Remove	Salected Attributes Sample Name Dot (,) Analysis Protocol Name Dot (,) Unique Time Stamp Integer

Figure 19 File name convention example

Figure 20 on page 164 shows the folders and files generated by the results group, file name convention, run name, and injections shown in Figure 17 on page 163, Figure 18 on page 163, and Figure 19 on page 164.

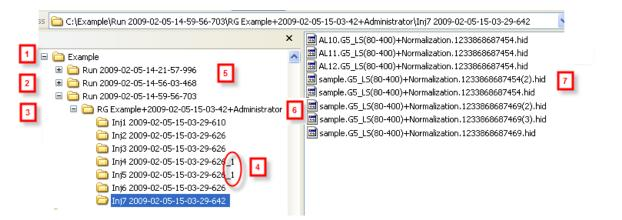


Figure 20 Folder hierarchy and file naming example

1	File location from results group <ul> <li>Custom file location</li> </ul>
2	Instrument Run Name folder from results group 🔽 Include an Instrument Run Name folder
3	Results group Name folder from results group 🔽 Include a Result Group Name folder
4	Injection folder from results group 🔽 Include an Injection folder
	Duplicate injections indicated with _n where n is the number of duplicates.
5	Run name (default or user-defined) from injection list Run Name: Run 2009-02-05-14-59-56-70:
6	Results group name syntax from results group
	RG Example+ <start date="" instrument="" run="" stamp="" time="">+<logged in="" name="" user=""></logged></start>
7	File name syntax from file name convention
	<sample name="">.<analysis name="" protocol="">.<unique integer="" stamp="" time=""></unique></analysis></sample>

# Instrument protocol library

#### Instrument protocol overview

An instrument protocol contains the parameters that control the instrument during data acquisition. An instrument protocol is a required element of an assay for all applications.

When you create an assay, you add one or more instrument protocols to the assay. If you add these items from the library, a *copy* of the items is added to the assay, and can be modified independently from the original items stored in the library. For information on how changes are tracked if auditing is enabled, see "Audit action" on page 210.

#### Create a new instrument protocol

If factory-provided instrument protocols do not suit your needs, you can create new instrument protocols:

- 1. Access the Instrument Protocols library.
- 2. Click 式 Create.
- 3. In the Create New Instrument Protocol dialog box (Figure 21 on page 166), select an application type: Sequencing, Fragment, or HID. The run module selection list is filtered based on the application you select.
- 4. Specify settings (Table 10 on page 166).
- 5. Save the assay:
  - If you are creating the assay from the Library, click **Save**.
- 블 3500 Data Collection Software Edit 🔻 Dashboard Library Maintenand 🗳 Library Resources 🖬 🖉 🖬 Filter: All 💥 Manage Plates Instrument Prot Assays 1 FragmentAnaly File Name Conventions 2 AR HID36 PC Results Group 3 AR HID36\_PO 🚹 Analyze 4 🛝 HID36\_PO 5 AR HID36 PO 6 AR HID36\_PO Dye Sets
- If you are creating the assay from the Assign Plate Contents screen, click **Apply to Plate** or **Save to Library**.

🔚 Create New Instrument Protocol
Setup an Instrument Protocol       Image: Setup an Instrument Protocol         Protocol Name "FragmentAnalysis50_POP7xl_1" already exists in the Library.       Image: Setup an Instrument Protocol
0
Application Type: Fragment 🔽 Capillary Length: 50 🔽 cm Polymer: POP7 🔽
Dye Set: G5
Instrument Protocol Properties
* Run Module: FragmentAnalysis50_POP7xI
* Protocol Name: FragmentAnalysis50_POP7xl_1
Description:
Oven Temperature (°C):       60       Run Voltage (kVolts):       19.5       PreRun Voltage (kVolts):       15       Injection Voltage (kVolts):       1.6         Run Time (sec.):       1330       PreRun Time (sec.):       180       Injection Time (sec.):       15       Data Delay (sec.):       1
▼ Advanced Options
Following values are not recommended to be changed.
Voltage Tolerance (kVolts):       0.7         Voltage # of Steps (nk):       30         Voltage Step Interval (sec.):       15
First Read Out Time (ms): 200 Second Read Out Time (ms): 200
Normalization Target: 4500.0 Normalization Factor Threshold Min: 0.3 Normalization Factor Threshold Max: 3.0
Close

Figure 21 Create New Instrument Protocol – normalization parameters circled in red are displayed for fragment analysis and HID applications only

# Instrument protocol settings

#### Table 10 Instrument protocol settings

Setting	Description
Application Type	Sequencing
	Fragment analysis
	• HID
Capillary Length, Polymer, Dye set	Capillary length, polymer type, and dye set with which the protocol will be used
Run module	Factory-provided modules that specify instrument control parameters. For more information, see "Run modules" on page 263.
Protocol name	Name of the protocol. Names must be unique.
Locked	When enabled, allows the entry to be unlocked and modified only by the user who created it, the administrator, or another user with unlock permissions. Useful when your system includes the SAE module (described in Chapter 7, "Use Security, Audit, and E-Sig Functions (SAE Module)" on page 197.
Description	Optional text entry.

Setting	Description
Oven temperature ( °C)	Temperature setting for main oven throughout run.
Run voltage (kVolts)	Final sample electrophoresis separation run voltage.
Prerun voltage (kVolts)	Pre run voltage setting before sample injection.
Injection voltage (kVolts)	Injection voltage setting for sample injection.
Run time (sec)	Length of time data is collected after voltage is ramped up to the run voltage and the run starts.
PreRun time (sec)	Prerun voltage time.
Injection time (sec)	Sample injection time.
Data delay (sec)	Time from the start of separation to the start of sample data collection.
Advanced options – Do not	change unless advised otherwise by Applied Biosystems support personnel
Voltage tolerance (kVolts)	Maximum allowed voltage variation.
Voltage # of Steps (nk)	Number of voltage ramp steps to reach Run Voltage.
Voltage step interval (sec)	Dwell time at each voltage ramp step.
First read out time (ms)	The interval of time for a data point to be produced. First ReadOut time should be equal to Second ReadOut time.
Second read out time (ms)	The interval of time for a data point to be produced. Second ReadOut time should be equal to First ReadOut time.
	<b>s only</b> : Normalization parameters – Leave at default settings (for information on how these eview normalized data" on page 90).
Normalization Target	The expected average RFU for the subset of peaks in the GS600 LIZ <sup>®</sup> v2 size standard used for normalization.
	The default value for each run module has been experimentally determined based on the average peak height of selected peaks in the GS600 size standard with a specific injection time.
	<b>IMPORTANT!</b> If you change the injection time in an instrument protocol, adjust the Normalization Target proportionately. For example, for an instrument protocol with an injection time of 10 seconds and a Normalization Target of 2000: if you change the injection time to 15 seconds (50% increase), change the Normalization Target to 3000 (50% increase).
Normalization Factor	The passing range for Normalization Factor (default range is 0.3 to 3.0).
Thresholds	<b>IMPORTANT!</b> Increasing the factor threshold above 3.0 may cause amplification of noise.
	If the calculated Normalization Factor is outside the Normalization Factor range, the software multiplies the peak heights of the sample by the low or high Normalization Factor threshold setting (for example, if the Normalization Factor range is 0.3 to 3.0 and the calculated Normalization Factor is 5, the software applies a Normalization Factor of 3.0).
	Average peak height of the subset of peaks in the GS600 LIZ® v2 size standard used
Normalization Factor	for normalization divided by the Normalization Target.

#### Table 10 Instrument protocol settings (continued)

# Dye sets library

### Dye set overview

A dye set defines the following for an instrument protocol:

- Dye color(s)
- Order of the dye peaks in the standard
- Spectral analysis parameters

When you create an instrument protocol, you add a dye set to the protocol. If you add this item from the library, a *copy* of the item is added to the assay, and can be modified independently from the original item stored in the library. For information on how changes are tracked if auditing is enabled, see "Audit action" on page 210.

### Create a new dye set

If factory-provided dye sets do not suit your needs, you can create new dye sets:

- 1. Access the Dye Sets library.
- 2. Click 📰 Create.
- 3. In the Create New Dye Set dialog box (Figure 22 on page 169).
- 4. Specify settings (Table 11 on page 169).
- 5. Click Save.



🔚 Create New Dye Set	×
Setup a Dye Set	Un.
🔇 Dye Set Name is a required field. Provide a unique value.	
	4
	0
	ocked
	ockea
* Chemistry Matrix Standard	
* Dye Set Template GS Template	
Arrange Dyes	
Calibration Peak Order 5 4 3 2 1	
▼ Parameters	
The parameters will be used for instruments configured with 50cm capillary array and polymer POP7	
Matrix Condition Number Upper Limit 13.5	
Locate Start Point * After Scan 1000 * Before Scan 5000	
* Limit Scans To 3250	
Sensitivity 0.4	
* Minimum Quality Score 0.95	
Notes	
Applied Biosystems	<b>A</b>
	<b>v</b>
Close	Save

Figure 22 Create New Dye Set

#### Table 11 Dye set settings

Setting	Description	
Dye Set Name	Name of the dye set. Names must be unique.	
Locked	When enabled, allows the entry to be unlocked and modified only by the user who created it, the administrator, or another user with unlock permissions. Useful when your system includes the SAE module (described in Chapter 7, "Use Security, Audit, and E-Sig Functions (SAE Module)" on page 197.	
Chemistry	The standard for which you are creating the dye set: Sequencing Standard or Matrix standard	
Dye Set Template	Factory-provided template upon which to base the dye set. The Any Dye template can be used for applications that do not use all of the dye colors contained in the matrix standard kits used for spectral calibration.	

#### Table 11 Dye set settings

Setting	Description
Arrange Dyes	Displays the dyes and the peak order for the dye set template selected.
	Editable only for AnyDye template:
	• Dye Selection – Specifies the dyes to use for calibration
	• Reduced Selection – Specifies the dyes used in the samples.
	For example, if you use the 5 dye kit and have samples with only blue peaks, you can "reduce" or deconvolute with blue and orange (size standard) dyes only.
Parameters	Specifies the Quality Value, Condition Number, Scan, and Sensitivity requirements for the dye set.
Notes	Optional text entry.

# Size standards library

#### Size standard overview

A size standard defines the sizes of known fragments. It is used to generate a standard curve. The standard curve is used to determine the sizing of unknown samples.

When you create a sizecalling (fragment) or QC (HID) protocol, you add a size standard to the protocol. If you add this item from the library, a *copy* of the item is added to the protocol, and can be modified independently from the original items stored in the library. For information on how changes are tracked if auditing is enabled, see "Audit action" on page 210.

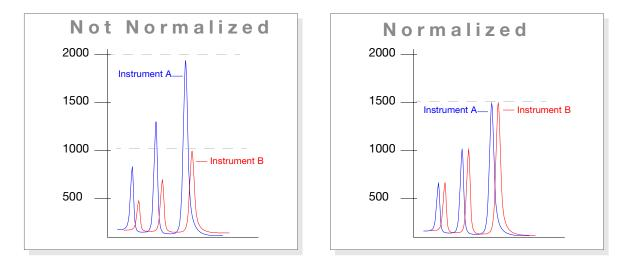
#### Normalization size standards provided

The library contains factory-provided normalized size standards that you can use to normalize fragment analysis and HID data:

- Fragment analysis:
  - GS600LIZ+Normalization
  - GS600(60-600)LIZ+Normalization For applications that have primer peaks that obscure the 20 and 40-mer peaks of the GS600 size standard.
- HID:
  - GS600(80-400)LIZ+Normalization

Normalization corrects for instrument, capillary, and injection variability. For each sample, the software calculates a normalization factor based on a threshold setting. The normalization factor is used as a multiplier to adjust the peak height of the sample peaks relative to the GS600  $LIZ^{(R)}$  v2.0 size standard peaks.

**IMPORTANT!** Normalization is not applied to samples with failing sizing quality. Select a size standard definition file appropriate for your application that accurately sizes samples. For example, if your application includes small fragments that may be obscured by primer peaks, or large fragments that may not be present due to slower migration rates, specify a size standard definition file that eliminates these fragments from sizing.

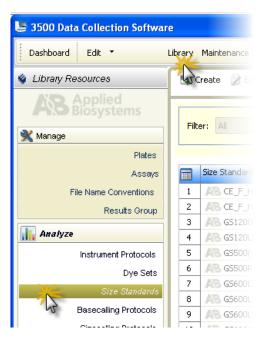


For more information, see "Review normalized data" on page 90.

### Create a new size standard

If factory-provided size standards do not suit your needs, you can create new size standards:

- 1. Access the Size Standards library.
- 2. Click 📑 Create.
- 3. In the Create New Size Standard dialog box (Figure 22 on page 169), enter a size standard name.
- 4. (Optional):
  - Select the Locked check box. When enabled, allows the entry to be unlocked and modified only by the user who created it, the administrator, or another user with unlock permissions. Useful when your system includes the SAE module (described in Chapter 7, "Use Security, Audit, and E-Sig Functions (SAE Module)" on page 197.



- Enter a description.
- 5. Select a dye color.
- 6. Enter sizes in the list on the left. Separate sizes with a comma, space, or return.
- 7. Click Add Sizes.

8. Click Save.

🔚 Create New Size Standard	×
Setup a Size Standard ⑧ Size Standard Name is a required field. Provide a unique value.	
* Size Standard Locked Description: * Dye Color:	Ø
Enter sizes in the field below separated by a comma, space, or return then click the "Add Size(s)>>" button to add them to the current size standard definition. Enter new Size Standard definition: (e.g. 11.0, 34.2, 55) Add Size(s) >>	Delete Selected Sizes
Close	Save

Figure 23 Create New Size Standard

# Modify a factory-provided normalization size standard

- 1. Select a factory-provided normalization size standard (indicated in the name with "+Normalization.")
- 2. Click 🕎 Duplicate.
- 3. Edit the copy of the normalized size standard. The size standard peaks used to normalize the data are displayed in gray and are not editable.
- 4. Click Save.

# Basecalling protocols library (primary analysis – sequencing)

#### Basecalling protocol overview

A basecalling protocol is the required primary analysis protocol for sequencing applications.

A basecalling protocol defines the settings used by the sequencing basecallers to assign base calls to each detected peak and assign a quality value:

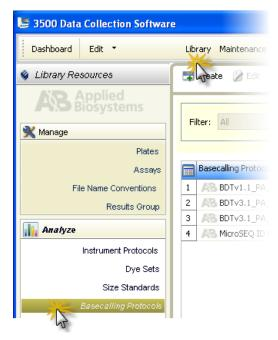
- Analysis settings
- Ranges for the sequencing quality flags displayed in View Results

When you create a sequencing assay, you add a basecalling protocol to the assay. If you add this item from the library, a *copy* of the item is added to the assay, and can be modified independently from the original item stored in the library. For information on how changes are tracked if auditing is enabled, see "Audit action" on page 210.

#### Create a new basecalling protocol

If factory-provided basecalling protocols do not suit your needs, you can create new basecalling protocols:

- 1. Access the Basecalling Protocols library.
- 2. Click 🛒 Create.
- In the Analysis Settings tab of the Create New Basecalling Protocol dialog box (Figure 24 on page 175), specify settings (see Table 12 on page 175).
- 4. Click **QV Settings**. In the QV Settings tab of the Create New Basecalling Protocol dialog box (Figure 25 on page 177), then specify settings and Table 13 on page 178).
- 5. Click Save.



堤 Create New Basecalling Protocol 🛛 🛛 🔀
Setup a Basecalling Protocol       Image: Constraint of the set of the
* Protocol Name:     Locked       Description:        Basecaller:     KB 1.4.1
Analysis Settings QV Settings
Mobility File     Quality Threshold     Mixed Bases Threshold     Mixed Bases Threshold     Analyzed Data Scaling     Clear Range Methods     Summary of current settings     Mobility File: KB_3500_POP7_BDTv3     Quality Threshold: Do not assign N's to Basecall     Mixed Base Threshold: 25.0%     Scaling: True Profile     Clear Range Methods: Use quality values.
Close Save

Figure 24 Create New Basecalling Protocol – Analysis Settings

Table 12	Basecalling protocol – Analysis settings
----------	--

Setting	Description		
Name	Name of the protocol. Names must be unique.		
Locked	When enabled, allows the entry to be unlocked and modified only by the user who created it, the administrator, or another user with unlock permissions. Useful when your system includes the SAE module (described in Chapter 7, "Use Security, Audit, and E-Sig Functions (SAE Module)" on page 197.		
Description	Optional text entry.		
Basecaller	Basecalling algorithm used to identify bases.		
	<b>Note:</b> The basecaller version listed in the basecalling protocol is a 3-digit number. The version listed in sequencing results is a 4-digit number. The fourth digit is an internal number used by the software.		
Mobility file	Compensates for mobility differences between dyes and primers, correcting the color code to the chemistry used to label the DNA during instrument processing.		

Table 12	Basecalling protocol - Anal	ysis settings (continued)
----------	-----------------------------	---------------------------

Setting	Description
Quality Threshold	Basecall Assignment (ambiguous bases):
	<ul> <li>Do not assign N's to basecalls</li> </ul>
	<ul> <li>Assign N's to basecalls with QV&lt;15 – Bases with a QV less than the threshold display N instead of the base letter</li> </ul>
	<ul> <li>Ending base – Last base on which to perform basecalling:</li> </ul>
	– At PCR Stop
	<ul> <li>After X number of Bases</li> </ul>
	<ul> <li>After X number of Ns in X number of Bases</li> </ul>
	<ul> <li>After X number of Ns</li> </ul>
	Note: If you have short PCR products, select the At PCR Stop check box.
Mixed bases threshold	When enabled, allows the software to determine the secondary peak height where the base position is considered a potential mixed base.
	Adjust this parameter by dragging the bar in the display or typing in a numeric value.
	Use Mixed Base Identification
	Do not assign a mixed base when the secondary peak height is $\leq to 25\%$
Analyzed Data Scaling	Determines scaling of the processed traces. This parameter does not affect the accuracy of the basecalling.
	• <b>True Profile</b> – The processed traces are scaled uniformly so that the average height of peaks in the region of strongest signal is about equal to a fixed value. The profile of the processed traces will be very similar to that of the raw traces.
	• Flat Profile – The processed traces are scaled semi-locally so that the average height of peaks in any region is about equal to a fixed value. The profile of the processed traces will be flat on an intermediate scale (> about 40 bases).
Clear range methods	• Use clear range minimum and maximum – Specifies the first and last base in the range to consider, or trims the specified number of bases from the 3' end.
	• Use quality values – Sets a window with a specified number of allowed low-quality bases by removing bases until there are < X number of bases per Z number of bases with QV < Y.
	<ul> <li>Use identification of N cells – Sets a window with a specified number of allowed ambiguous base calls (Ns) by removing bases until there are &lt; X number of Ns per Y number of bases.</li> </ul>

블 Create New Basecalli	ng Protocol			×
Setup a Basecalling Pro		unique value.		
Protocol Name: Description: Basecaller: KB 1.4.				
Analysis Settings QV Settin	ngs Fail If Value Is	Sucpert Dange	0	
Contiguous Read Length Trace Score QV20+	<ul> <li>100</li> <li>15</li> <li>100</li> </ul>	Suspect Range 100-300 15-30 100-300	Pass if Value Is           >=         300           >=         30           >=         300	
Close	(	Save		

Figure 25 Create New Basecalling Protocol – QV Settings

QV settings are quality value ranges used in the following screens:

- Monitor Run screen The state of the QV flag:
  - If all three values are in the pass range, the QV flag in Monitor Run is set to
     (green).
  - If any values are in the suspect range, the QV flag in Monitor Run is set to
     (yellow).
  - If any value fails are in the fail range, the QV flag in Monitor Run is set to
     (red).
- View Sequencing Results > Metric Analysis Results table The pass/check/fail status for Trace Score Quality, CRL Quality, and QV20+ Quality results.

#### Table 13 Basecalling protocol – QV settings

Setting	Description
Contiguous Read Length	The longest uninterrupted segment of bases with an average Quality Value (QV) $\ge$ 20.
	In addition to evaluating the QV of a base call, the software considers the QV of adjacent bases within a $\pm$ 20-bp moving average to determine a contiguous read length based on quality values: the software starts from the 5' end and calculates the average QV across a moving window size of 20, sliding 1 bp at a time, to the 3' end. The resulting longest contiguous segment is determined as the CRL.
Trace Score	The average basecall quality value (QV) of bases in the clear range sequence of a trace.
QV20+	The total number of bases in the entire trace with quality values $\geq 20$ .

# Sizecalling protocols library (primary analysis - fragment)

### Sizecalling protocol overview

A sizecalling protocol is the required primary analysis protocol for fragment applications.

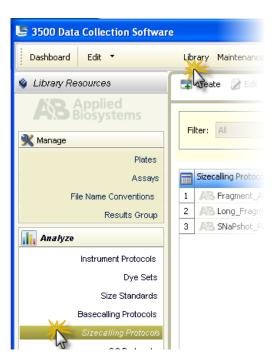
A sizecalling protocol defines peak detection, sizing, and quality values.

When you create a fragment assay, you add a sizecalling protocol to the assay. If you add this item from the library, a *copy* of the item is added to the assay, and can be modified independently from the original item stored in the library. For information on how changes are tracked if auditing is enabled, see "Audit action" on page 210.

# Create a new sizecalling protocol

If factory-provided sizecalling protocols do not suit your needs, you can create new sizecalling protocols:

- 1. Access the Sizecalling Protocols library.
- 2. Click 📑 Create.
- In the Analysis Settings tab of the Create New Sizecalling Protocol dialog box (Figure 26 on page 180), specify settings (see Table 14 on page 180).
- 4. Click **QC Settings**. In the QC Settings tab of the Create New Sizecalling Protocol dialog box (Figure 27 on page 183), then specify settings and Table 15 on page 183).
- 5. Click Save.



Create New S	izecalling Protocol				
Setup a Sizecal Setup of Sizecal	ling Protocol a required field. Provide	e a unique value.			<b>K</b>
* Protocol Name:					Locked
Description:					
Size Standard:	G5600LIZ	*			
Sizecaller:	SizeCaller v1.1.0 🔽				
Analysis Settings	QC Settings				
					0
Analysis Range:	Full 💌	Sizing Range: Full	×	Size Calling Method:	Local Southern 🔽
Analysis Start Po		Sizing Start Size: 0	- Hulloan	Primer Peak:	Present 🔽
Analysis Stop Po		Sizing Stop Size: 100	000		
	<b>∣</b> Blue	Green Vell	ow 🔽 Red	V Purple	🗸 Orange
Minimum Da alu					
Minimum Peak	Height 175	175 175	5 175	175	175
Common Settin	ngs				
		Use Smoothing	None 😽		
	Use Bas	elining (Baseline Window (Pts))	51		
		Minimum Peak Half Width	2		
		Peak Window Size	15		
		Polynomial Degree	3		
		51 55			
		Slope Threshold Peak Start	0.0		
		Slope Threshold Peak End	0.0		
					]
Close					Save

Figure 26 Create New Sizecalling Protocol – Analysis Settings

**IMPORTANT!** Normalization is not applied to samples with Size Quality flags. Specify analysis settings that accurately detect and size the size standard, and QC settings with appropriate pass fail ranges. The 3500 Series Data Collection Software does not support re-analyzing data with new settings.

#### Table 14 Sizecalling protocol – Analysis settings

Setting	Description		
Protocol Name	Name of the protocol. Names must be unique.		
Description	Optional text entry.		
Size standard	Size standard definition in the software that corresponds to the dye set used in the chemistry.		
	To apply normalization, select a normalization size standard (see "Normalization size standards provided" on page 171).		

Setting	Description
Analysis Range	Specify the range (in data points) to analyze:
	• <b>Full Range</b> to analyze the entire scan region as collected by the genetic analysis instrument, including the primer peak.
	• <b>Partial Range</b> to analyze only data points within a specified range. Enter Start Point in data points after the primer peak and before the first required size standard peak Enter a Stop Point after the last required size standard fragment. Start and Stop points may vary from instrument to instrument and platform to platform. Display raw data to determine the appropriate analysis range.
	Data points outside the specified analysis range are ignored.
	<b>Note:</b> Ensure the Analysis Range contains all size standard fragments included in the Sizing Range specified below.
Sizing Range	Specify the size range (in base pairs) appropriate for the kit you are using:
	• All Sizes for the software to analyze fragments of all sizes in the Analysis Range.
	• Partial Sizes for the software to analyze only fragments within a specified range. Enter a Start Size and a Stop Size appropriate for the size standard used.
Size Calling Method	• Local Southern - (default) Determines the fragment sizes using the reciprocal relationship between fragment length and electrophoretic mobility.
	3rd Order Least Squares - Uses regression analysis to build a best-fit size calling curve.
	<ul> <li>2nd Order Least Squares - Uses regression analysis to build a best-fit size calling curve.</li> </ul>
	<ul> <li>Cubic Spline Interpolation - Forces the sizing curve through all the known points of the selected size standard.</li> </ul>
	Global Southern Method - Compensates for standard fragments with anomalous electrophoretic mobility (similar to least squares methods).
Primer Peak	If the primer peaks in your application obscure peaks of interest, select <b>Present</b> . Selecting Present instructs the algorithm to ignore primer peaks. Primer peaks are stil displayed in the trace.
	<b>Note:</b> If this setting does not allow detection of the 20 and 40-mer peaks for samples that use the GS600 LIZ size standard, running samples with the GS600(60-600)LIZ+Normalization may allow detection of the peaks.
Peak Amplitude Thresholds	Specify the threshold (RFU) for peak detection for each dye color. Peaks below the threshold are not detected.
	For example, if you use the default values of 175, peaks with heights equal to or greater than 175 are detected. Peaks with heights below 175 are still displayed in the electropherogram plots but are not detected or labeled.
	<b>Note:</b> Set the peak amplitude thresholds to 175 in your GeneMapper <sup>®</sup> Software analysis method.
Smoothing	Select an option to smooth the outline of peaks and reduce the number of false peaks detected:
	• <b>None</b> (default) to apply no smoothing. Best if the data display sharp, narrow peaks of interest.
	• Light to provide the best results for typical data. Light smoothing slightly reduces peak height.
	<ul> <li>Heavy for data with very sharp, narrow peaks of interest. Heavy smoothing can significantly reduce peak height.</li> </ul>

#### Table 14 Sizecalling protocol – Analysis settings (continued)

#### Table 14 Sizecalling protocol – Analysis settings (continued)

Setting	Description
Baseline Window	Specify a window to adjust the baseline signals of all detected dye colors to the same level for an improved comparison of relative signal intensity. Note the following:
	<ul> <li>A small baseline window relative to the width of a cluster, or grouping of peaks spatially close to each other, can result in shorter peak heights.</li> </ul>
	<ul> <li>Larger baseline windows relative to the peaks being detected can create an elevated baseline, resulting in peaks that are elevated or not resolved to the baseline.</li> </ul>
Min. Peak Half Width	Specify the smallest half peak width at full height for peak detection. The range is 2 to 99 data points.
Polynomial Degree	Polynomial Degree cannot be greater than Peak Window Size.
	Adjust to affect the sensitivity of peak detection. You can adjust this parameter to detect a single base pair difference while minimizing the detection of shoulder effects and/or noise.
	The peak detector calculates the first derivative of a polynomial curve fitted to the data within a window that is centered on each data point in the analysis range.
	Using curves with larger polynomial degree values allows the curve to more closely approximate the signal and, therefore, captures more of the peak structure in the electropherogram.
	For information on optimizing Polynomial Degree and Peak Window Size, see
Peak Window Size	Enter a window width in data points for peak detection sensitivity. If more than one peak apex is within the window, all are labeled as a single peak. Note the following:
	The maximum value is the number of data points between peaks.
	The Peak Window Size setting is limited to odd numbers.
	To increase peak detection sensitivity: Increase polynomial degree, decrease peak window size.
	To decrease peak detection sensitivity: Decrease polynomial degree, increase peak window size.
Slope Thresholds Peak Start and End	• <b>Peak Start</b> - The peak starts when the first derivative (slope of the tangent) in the beginning of the peak signal before the inflection point becomes equal to or exceeds the "Peak Start" value. This threshold is set to 0 by default, which means that the peak will normally start at the leftmost point where the slope of the tangent is closest to 0° (horizontal line). A value other than 0 moves the peak start point toward its center. The value entered must be non-negative.
	• <b>Peak End</b> - The peak ends when the first derivative (slope of the tangent) in the end of the peak signal after the inflection point becomes equal to or exceeds the "Peak End" value. This value is set to 0 by default, which means that the peak will normally end at the rightmost point where the slope of the tangent is closest to 0° (horizontal line). A value other than 0 moves the peak end point toward its center. The value entered in this field must be non-positive.

Protocol Name is a required f		Let a
nalysis Settings QC Settings	5	
		G
Size Quality	272	
Fail if Value is	Suspect Range	Pass if Value is
< 0.25	0.25 - 0.75	≥ 0.75
Assume Linearity from (bp):	0 To (bp): 80	0
Pull Up		
Actuate Pull-Up flag if Pull-Up	Ratio ≤ 0.05 and Pul	l-Up Scan ≤ 1

Figure 27 Sizecalling Protocol – QC Settings

**IMPORTANT!** Normalization is not applied to samples with Size Quality flags. The 3500 Series Data Collection Software does not support re-analyzing data with new settings.

Table 15 Sizecalling protocol – QC settings

Setting	Description
Size Quality	Enter the Pass Range and the Low Quality Range for the SQ flag displayed in View Fragment Results.
	Results that are within the Pass range are flagged as [] (Pass). Results that are within the Low Quality range are flagged as () (Low Quality). Results that are between the Pass and Low Quality ranges are flagged (Check).
	For example, with a Pass Range of 0.75 to 1.0 and a Low Quality Range of 0.0 to 0.25, any result above 0.75 is, any result at 0.25 or lower is, and any result between 0.26 to 0.74 is
	How Size Quality is determined
	The Size Quality algorithm evaluates the similarity between the fragment pattern for the size standard dye specified in the size standard definition and the actual distribution of size standard peaks in the sample, calculates an interim SQ (a value between 0 and 1).
Assume Linearity	Defines the expected linear range. Useful in large fragment size standards where non- linearity might be expected.
Pull-Up	Enter the pull-up ratio and tolerance for pull-up peak identification.
	A pull-up peak is identified when the peak height of the minor peak is:
	• $\leq$ X% (pull-up ratio) of the major peak <i>and</i>
	• Within ±Y data point (pull-up scan) of the major peak
	When at least one peak is identified as a pull-up peak, the 📥 (Check) flag is displayed for the Spectral Pull-Up quality flag in View Fragment Results.

# QC protocols library (primary analysis - HID)

### QC protocol overview

A QC protocol is the required primary analysis protocol for HID applications. A QC protocol defines peak detection, sizing, and quality values.

When you create an HID assay, you add a QC protocol to the assay. If you add this item from the library, a *copy* of the item is added to the assay, and can be modified independently from the original item stored in the library. For information on how changes are tracked if auditing is enabled, see "Audit action" on page 210.

### Create a new QC protocol

If factory-provided QC protocols do not suit your needs, you can create new QC protocols:

1. Access the QC Protocols library.

블 3500 Data Collection Softwar	e
Dashboard Edit 🔻	Library Maintenance Tools
🗳 Library Resources	📑 🤯ate 🖉 Edit 🖽 Dur
AB Applied Biosystems	Filter: All
Plates Assays	QC Protocol Name
File Name Conventions	1 AB F_LS(75-400)
Results Group	2 🗛 F_LS(75-450)
	3 AB G5_3rd(80-400)+Nor
Analyze	4 AB G5_3rd(80-400)
Instrument Protocols	5 AB G5_LS(80-400)+Nor.
Dye Sets	6 🕂 G5_LS(80-400)
Size Standards	
Basecalling Protocols	
Sizecalling Protocols	
QC Protocols	
Saguande Analysis Drotocols	1. I.

- 2. Click 式 Create.
- 3. In the Analysis Settings tab of the Create New QC Protocol dialog box (Figure 28 on page 185), specify settings (see Table 16 on page 185).
- 4. Click **QC Settings**. In the QC Settings tab of the Create New QC Protocol dialog box (Figure 29 on page 188), specify settings (Table 17 on page 188).
- 5. Click Save.

Create New C	QC Prote	ocol					
etup a QC Pro Protocol Name i		ed field. Provid	e a unique value.				
* Protocol Name:	Î						
escription:							
iize Standard:	G5600_	_LIZ+Normaliza	tion_(80-400) 🔽				
iizecaller:	SizeCal	ler v1.1.0 🔽	]				
Analysis Settings	QC Set	tings					
Analysis Range: Analysis Start P Analysis Stop Po	oint: 0	Number of	Sizing Range Sizing Start S Sizing Stop S	iize: 0		Size Calling Method:	Local Southern 💊
Peak Ar	mplitude	<b>Blue</b> 175	<b>Green</b>	Yellow	<b>Red</b>	Purple	Orange
Common Setti							
		Use Baseli	Use Si ning (Baseline Windo	moothing Light			
			Minimum Peak H	alf Width			
			Peak Win	dow Size	5		
			Polynomia	al Degree			
			Slope Threshold P		.0		
			Slope Threshold I		.0		
<u></u>							

Figure 28 Create New QC Protocol – Analysis Settings

**IMPORTANT!** The default values in the QC protocol templates (other than peak amplitude threshold values) have been optimized for each kit. You must optimize and validate peak amplitude threshold values during internal HID validation. If you modify other settings, ensure that the size standard is accurately detected and sized with the new settings. Normalization is not applied to samples with  $\mathbf{X}$  Size Quality flags. The 3500 Series Data Collection Software does not support re-analyzing data with new settings.

Table 16	QC protocol	<ul> <li>Analysis</li> </ul>	settings
----------	-------------	------------------------------	----------

Setting	Description
Protocol Name	Name of the protocol. Names must be unique.
Description	Optional text entry.

Table 16 Q	C protocol – Analy	sis settings	(continued)
------------	--------------------	--------------	-------------

Setting	Description
Size standard	Size standard definition in the software that corresponds to the dye set used in the chemistry.
	To apply normalization, select a normalization size standard (see "Normalization size standards provided" on page 171).
Analysis Range	Select <b>Full</b> to collect data points for the entire scan region, including the primer peak. You can specify a limited analysis range in the GeneMapper <sup>®</sup> <i>ID-X</i> Software.
	Note: If you select Partial, ensure that the Analysis Range contains all size standard fragments included in the Sizing Range specified below.
Sizing Range	Select <b>Partial</b> , then specify <b>80</b> to <b>400</b> to limit the fragment sizes evaluated for the size standard.
	If you specify sizes outside this range, the Sizing Quality may fail.
Size Calling Method	Select the method to determine the molecular length of unknown fragments appropriate for the AmpF/STR <sup>®</sup> kit you use:
	<ul> <li>Local Southern - (default) Determines the fragment sizes using the reciprocal relationship between fragment length and electrophoretic mobility. The unknown fragment is surrounded by two known-sized fragments above and one below, then two below and one above. The results are averaged and the size of the allele is determined</li> </ul>
	<ul> <li>Identifiler<sup>®</sup> kit</li> <li>SEfiler Plus<sup>™</sup> kit</li> <li>Sinofiler<sup>™</sup> kit</li> <li>Yfiler<sup>®</sup> kit</li> <li>SGM Plus<sup>®</sup> kit</li> </ul>
	<ul> <li>3rd Order Least Squares - Uses regression analysis to build a best-fit size calling curve</li> <li>MiniFiler<sup>™</sup> reagent.</li> </ul>
	Size calling options for kits other than those listed above are:
	<ul> <li>2nd Order Least Squares - Uses regression analysis to build a best-fit size calling curve.</li> </ul>
	<ul> <li>Cubic Spline Interpolation - Forces the sizing curve through all the known points of the selected size standard.</li> </ul>
	• <b>Global Southern Method</b> - Compensates for standard fragments with anomalous electrophoretic mobility (similar to least squares methods).
with the new settings. N Collection Software doe	odify peak detection settings, ensure that the size standard is accurately detected and sized lormalization is not applied to samples with Size Quality flags. The 3500 Series Data es not support re-analyzing data with new settings. For more information on peak detection neMapper <sup>®</sup> <i>ID-X</i> Software Reference Guide.
Smoothing	Select an option to smooth the outline of peaks and reduce the number of false peaks detected:
	• None to apply no smoothing. Best if the data display sharp, narrow peaks of interest.
	• Light (default) to provide the best results for typical data. Light smoothing slightly reduces peak height.
	• <b>Heavy</b> for data with very sharp, narrow peaks of interest. Heavy smoothing can significantly reduce peak height.
Baseline Window	Specify a window to adjust the baseline signals of all detected dye colors to the same leve for an improved comparison of relative signal intensity. Note the following:
	• A small baseline window relative to the width of a cluster, or grouping of peaks spatially

• Larger baseline windows relative to the peaks being detected can create an elevated baseline, resulting in peaks that are elevated or not resolved to the baseline.

Setting	Description
Peak Amplitude	<b>IMPORTANT!</b> Optimize these thresholds during internal HID validation.
Thresholds	Specify the threshold (RFU) for peak detection for each dye color. Peaks below the threshold are not detected.
	For example, if you use the default values of 175, peaks with heights equal to or greater than 175 are detected. Peaks with heights below 175 are still displayed in the electropherogram plots but are not detected or labeled.
	<b>Note:</b> Ensure that the same peak amplitude thresholds are used in secondary analysis software such as GeneMapper <sup>®</sup> (v4.1 or later)
Min. Peak Half Width	Specify the smallest half peak width at full height for peak detection. The range is 2 to 99 data points.
Polynomial Degree	Adjust to affect the sensitivity of peak detection. You can adjust this parameter to detect a single base pair difference while minimizing the detection of shoulder effects and/or noise
	The peak detector calculates the first derivative of a polynomial curve fitted to the data within a window that is centered on each data point in the analysis range.
	Using curves with larger polynomial degree values allows the curve to more closely approximate the signal and, therefore, captures more of the peak structure in the electropherogram.
Peak Window Size	Enter a window width in data points for peak detection sensitivity. If more than one peak apex is within the window, all are labeled as a single peak. Note the following:
	The maximum value is the number of data points between peaks.
	The Peak Window Size setting is limited to odd numbers.
	To increase peak detection sensitivity: Increase polynomial degree, decrease peak window size.
	To decrease peak detection sensitivity: Decrease polynomial degree, increase peak window size.
Slope Thresholds Peak	Not recommended for use with AmpFt/STR <sup>®</sup> kit data.
Start and End	• <b>Peak Start</b> - The peak starts when the first derivative (slope of the tangent) in the beginning of the peak signal before the inflection point becomes equal to or exceeds th "Peak Start" value. This threshold is set to 0 by default, which means that the peak will normally start at the leftmost point where the slope of the tangent is closest to 0° (horizontal line). A value other than 0 moves the peak start point toward its center. The value entered must be non-negative.
	<ul> <li>Peak End - The peak ends when the first derivative (slope of the tangent) in the end of the peak signal after the inflection point becomes equal to or exceeds the "Peak End" value. This value is set to 0 by default, which means that the peak will normally end at the rightmost point where the slope of the tangent is closest to 0° (horizontal line). A value other than 0 moves the peak end point toward its center. The value entered in th field must be non-positive.</li> </ul>

Table 16	QC protocol – Analysis settings (continued)
----------	---

Protocol Name:		
scription:		
e Standard:	G5600LIZ	
ecaller:	SizeCaller v1.1.0 💙	
alysis Settings	QC Settings	
		0
Size Quality Fail if Va	alue is Suspect Range Pass if Value is	

Figure 29 Create New QC Protocol – Analysis Settings

**IMPORTANT!** Normalization is not applied to samples with Size Quality flags. The 3500 Series Data Collection Software does not support re-analyzing data with new settings.

#### Table 17 QC Protocol – QC Settings

Setting	Description		
Size Quality	Enter the Pass Range and the Low Quality Range for the SQ flag displayed in View HID Results.		
	Results that are within the Pass range are flagged as [26] (Pass). Results that are within the Low Quality range are flagged as (26) (Low Quality). Results that are between the Pass and Low Quality ranges are flagged (26) (Check).		
	For example, with a Pass Range of 0.75 to 1.0 and a Low Quality Range of 0.0 to 0.25 any result above 0.75 is any result at 0.25 or lower is , and any result between 0.26 to 0.74 is		
Size Quality	How Size Quality is determined		
	The Size Quality algorithm evaluates the similarity between the fragment pattern for the size standard dye specified in the size standard definition and the actual distribution of size standard peaks in the sample, calculates an interim SQ (a value between 0 and 1)		
	Weighting		
	The Broad Peak (BD) threshold specified in the QC Protocol - QC Settings tab affects the SQ. To determine the final SQ value, the software:		
	• Evaluates size standard peak widths in the sample in the dye color specified in the size standard definition.		
	<ul> <li>If the width of any size standard peak in the sizing range exceeds the broad peak threshold, applies a 0.5 weighting factor: Interim SQ × (1–0.5)</li> </ul>		
	<b>Note:</b> The GeneMapper <sup>®</sup> <i>ID-X</i> Software allows you to set broad peak weighting. For more information, see the GeneMapper <sup>®</sup> <i>ID-X</i> Software Reference Guide.		
Broad Peak	Enter the maximum peak width (in base pairs).		
	When a peak width is greater than the threshold, the 🐣 (Check) flag is displayed for the BD (Broad Peak) quality flag in View HID Results.		

# Sequencing analysis protocols library (secondary analysis)

### Sequencing analysis protocol overview

A sequencing protocol is the optional secondary analysis (auto-analysis) protocol for SeqScape<sup>®</sup> Software v2.7 or later sequencing applications.

A sequencing analysis protocol defines the:

- Secondary analysis software (SeqScape<sup>®</sup> Software) location
- SeqScape<sup>®</sup> Software project, template, and specimen to use for auto-analysis

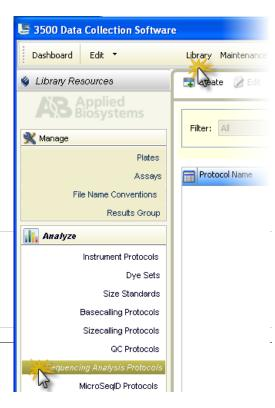
When you create a sequencing assay, you can optionally add a sequencing analysis protocol to the assay. If you add this item from the library, a *copy* of the item is added to the assay, and can be modified independently from the original item stored in the library. For information on how changes are tracked if auditing is enabled, see "Audit action" on page 210.

#### Create a new sequencing analysis protocol

- 1. Access the Sequencing Analysis Protocols library.
- 2. Click 式 Create.
- In the Create New Sequencing Analysis Protocol dialog box (Figure 30 on page 190), specify settings (see Table 18 on page 190).
- 4. Select the remaining secondary analysis items, then click **Save**.

**Note:** If the project, project template, or specimen of interest is not displayed in a list, re-select the secondary analysis software instance to update the list.

**IMPORTANT!** The auto-analysis settings you specify for the plate to run with this protocol must contain the same secondary



software and location settings. For more information, see "Create a new plate" on page 144.

etup a Sequencing Analysis Pr Protocol Name is a required field. Pro		
Protocol Name:		
Description:		
Application Type:	Sequencing	¥
* Secondary Analysis Software:	SeqScape	Ŧ
Secondary Analysis Sortware.		
Secondary Analysis Software Instance:	SeqScape_sanghahs-pc	•
	SeqScape_sanghahs-pc	•
Secondary Analysis Software Instance: Properties	SeqScape_sanghahs-pc	•
Secondary Analysis Software Instance: Properties	p53_v2	•

Figure 30 Create New Sequencing Analysis Protocol

Table 18	Create New Sequencing Analysis Protocol
----------	---

Setting	Description		
Protocol Name	Name of the protocol. Names must be unique.		
Description	Optional text entry.		
Lock	When enabled, allows the entry to be unlocked and modified only by the user who created it, the administrator, or another user with unlock permissions. Useful when your system includes the SAE module (described in Chapter 7, "Use Security, Audit, and E-Sig Functions (SAE Module)" on page 197.		
Application Type	Automatically set to Sequencing.		
Secondary Analysis Software	<b>IMPORTANT!</b> The secondary analysis software must be installed and properly configure with the 3500 Series Data Collection Software before it is listed as a selection in this screen. For information on setting up the SeqScape® Software for auto-analysis.		
Secondary Analysis Software Instance	Computer on which the secondary analysis software is running.		
Project	SeqScape software project to create.		
Project Template	Project template to use.		
Specimen	Specimen in which to save the sample data files.		
	Note: For each specimen a sequencing analysis protocol is required.		

# MicroSeq<sup>®</sup> ID protocols library (secondary analysis)

# MicroSeq<sup>®</sup> ID analysis protocol overview

A MicroSeq<sup>®</sup> ID protocol is the optional secondary analysis (auto-analysis) protocol for MicroSeq<sup>®</sup> ID Analysis Software v2.2 or later sequencing applications.

A MicroSeq<sup>®</sup> ID analysis protocol defines the:

- Secondary analysis software (MicroSeq® ID Analysis Software) location
- MicroSeq<sup>®</sup> ID Analysis Software project and specimen to use for auto analysis

When you create a sequencing assay, you can optionally add a MicroSeq<sup>®</sup> ID analysis protocol to the assay. If you add this item from the library, a *copy* of the item is added to the assay, and can be modified independently from the original item stored in the library. For information on how changes are tracked if auditing is enabled, see "Audit action" on page 210.

# Create a new MicroSeq® ID analysis protocol

- Access the MicroSeq<sup>®</sup> ID Protocols library.
- 2. Click 式 Create.
- In the Create New MicroSeq<sup>®</sup> ID Protocol dialog box (Figure 31 on page 192), specify settings (see Table 19 on page 192).
- 4. Select the remaining secondary analysis items, then click **Save**.

**Note:** If the project or specimen of interest is not displayed in a list, re-select the secondary analysis software instance to update the list.

**IMPORTANT!** The auto-analysis settings you specify for the plate to run with this protocol must contain the same secondary software and location settings. For more information, see "Create a new plate" on page 144.



블 Create New Mi	croSeqID Protocol			
Setup a MicroSe	Setup a MicroSeq® ID Protocol			
🔇 Protocol Name is a	a required field. Provid	e a unique value.		
* Protocol Name:				
Description:				
	Application Type:	Sequencing	~	
* Secondar	y Analysis Software:		~	
* Secondary Analysis	Software Instance:		~	
-Properties				
	* Project:		~	
	* Specimen:		~	
Close				

Figure 31 Create New MicroSeq<sup>®</sup> ID Protocol

Table 19	MicroSeq®	<b>ID</b> Analysis	protocol settings
----------	-----------	--------------------	-------------------

Setting	Description
Protocol Name	Name of the protocol. Names must be unique.
Description	Optional text entry.
Lock	When enabled, allows the entry to be unlocked and modified only by the user who created it, the administrator, or another user with unlock permissions. Useful when your system includes the SAE module (described in Chapter 7, "Use Security, Audit, and E-Sig Functions (SAE Module)" on page 197.
Application Type	Automatically set to Sequencing.
Secondary Analysis Software	<b>IMPORTANT!</b> The secondary analysis software must be installed and properly configured with the 3500 Series Data Collection Software before it is listed as a selection in this screen. For information on setting up the MicroSeq <sup>®</sup> <i>ID</i> Analysis Software for auto-analysis, see the <i>MicroSeq<sup>®</sup> ID</i> Analysis Software Getting Started <i>Guide</i> .
Secondary Analysis Software Instance	Computer on which the secondary analysis software is running.
Project	MicroSeq <sup>®</sup> ID software project and specimen to create.
Project Template	Project template to use.
Specimen	Specimen in which to save the sample data files.

# Fragment analysis protocols library (secondary analysis)

## Fragment analysis protocol overview

A fragment analysis protocol (GeneMapper<sup>®</sup> protocol) is the optional secondary analysis (auto-analysis) protocol for GeneMapper<sup>®</sup> Software v4.1 or later fragment applications.

A fragment analysis protocol defines the:

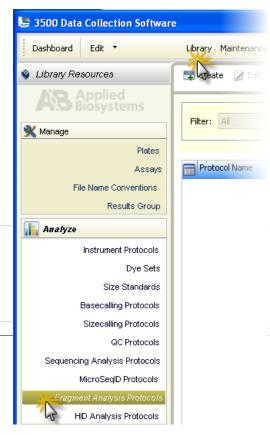
- Secondary analysis software (GeneMapper<sup>®</sup> Software) location
- GeneMapper<sup>®</sup> software analysis method, size standard, and panel that the GeneMapper<sup>®</sup> software will use during auto-analysis

## Create a new fragment analysis protocol

- 1. Access the Fragment Analysis Protocols library.
- 2. Click 📑 Create.
- In the Create New Fragment Analysis Protocol dialog box (Figure 32 on page 194), specify settings (see Table 20 on page 194).
- 4. Select the remaining secondary analysis items, then click **Save**.

**Note:** If the analysis method, size standard, or panel of interest is not displayed in a list, re-select the secondary analysis software instance to update the list.

**IMPORTANT!** The auto-analysis settings you specify for the plate to run with this protocol must contain the same secondary software and location settings. For more information, see "Create a new plate" on page 144.



Getup a GeneM	opper® Protocol		41
Protocol Name is	a required field. Provide a unique value.		
			0
* Protocol Name:			
Description:			
	Application Type: Fragment Analysis	~	
* Seconda	ry Analysis Software:	×	
* Secondary Analy:	is Software Instance:	~	
Properties			
	* Analysis Method:	*	
	* Size Standard:	~	
	Panel:	~	

#### Figure 32 Create New Fragment Analysis Protocol

Table 20	Fragment Analy	ysis protoco	settings
----------	----------------	--------------	----------

Setting	Description		
Protocol Name	Name of the protocol. Names must be unique.		
Description	Optional text entry.		
Lock	When enabled, allows the entry to be unlocked and modified only by the user who created it, the administrator, or another user with unlock permissions. Useful when your system includes the SAE module (described in Chapter 7, "Use Security, Audit, and E-Sig Functions (SAE Module)" on page 197.		
Application Type	Automatically set to Fragment analysis		
Secondary Analysis Software	<b>IMPORTANT!</b> The secondary analysis software must be installed and properl configured with the 3500 Series Data Collection Software before it is listed as selection in this screen.		
Secondary Analysis Software Instance	Computer on which the secondary analysis software is running		
Properties	GeneMapper^ $^{\ensuremath{\mathbb{R}}}$ software analysis method, size standard, and panel to use for autoanalysis.		

# HID analysis protocols library (secondary analysis)

## HID analysis protocol overview

An HID analysis protocol (GeneMapper<sup>®</sup> ID-X protocol) is the optional secondary analysis (auto-analysis) protocol for GeneMapper<sup>®</sup> ID-X Software v1.2 or later for HID applications.

An HID analysis protocol defines the:

- Secondary analysis software (GeneMapper<sup>®</sup> *ID-X* Software) location
- GeneMapper<sup>®</sup> *ID-X* Software analysis method, size standard, and panel that the GeneMapper<sup>®</sup> *ID-X* Software will use during auto-analysis

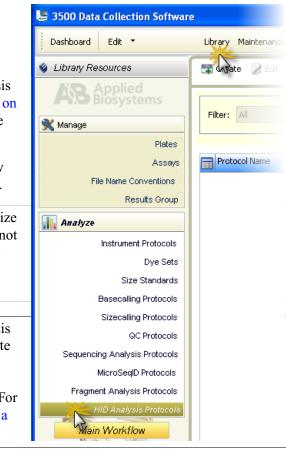
When you create an HID assay, you can optionally add an HID analysis protocol to the assay. If you add this item from the library, a *copy* of the item is added to the assay, and can be modified independently from the original item stored in the library. For information on how changes are tracked if auditing is enabled, see "Audit action" on page 210.

### Create a new HID analysis protocol

- 1. Access the HID Analysis Protocols library.
- 2. Click 📑 Create.
- 3. In the Create New HID Analysis Protocol dialog box (Figure 33 on page 196), specify settings (see Table 21 on page 196).
- 4. Select the remaining secondary analysis items, then click **Save**.

**Note:** If the analysis method, size standard, or panel of interest is not displayed in a list, re-select the secondary analysis software instance to update the list.

**IMPORTANT!** The auto-analysis settings you specify for the plate to run with this protocol must contain the same secondary software and location settings. For more information, see "Create a new plate" on page 144.



	apper® ID-X Protocol : a required field. Provide a unique value.	
		Y
		0
* Protocol Name:		
Description:		
	Application Type: HID	
* Seconda	ary Analysis Software:	
* Secondary Analy:	sis Software Instance:	
Properties		
	* Analysis Method:	
	* Size Standard:	
	Panel:	

Figure 33 Create New HID Analysis Protocol

Table 21	HID Anal	ysis protocol	l settings
----------	----------	---------------	------------

Setting	Description		
Protocol Name	Name of the protocol. Names must be unique.		
Description	Optional text entry.		
Lock	When enabled, allows the entry to be unlocked and modified only by the user who created it, the administrator, or another user with unlock permissions. Useful when your system includes the SAE module (described in Chapter 7, "Use Security, Audit, and E-Sig Functions (SAE Module)" on page 197.		
Application Type	Automatically set to HID.		
Secondary Analysis Software <b>IMPORTANT!</b> The secondary analysis software must be install configured with the 3500 Series Data Collection Software before selection in this screen. For information on setting up the Genel Software for auto-analysis, see the GeneMapper <sup>®</sup> <i>ID-X</i> Software <i>Guide</i> .			
Secondary Analysis Software Instance	Computer on which the secondary analysis software is running.		
Properties	GeneMapper <sup>®</sup> <i>ID-X</i> analysis method, size standard, and panel to use for auto-analysis.		

**Note:** If you are running a stand-alone version of the 3500 Series Data Collection Software (a version that is not installed on the instrument computer), you can create plates and protocols, then export them for use on the instrument computer.

# Section 1 Administrators

# Administrators overview of system security, auditing, and electronic signature

The SAE (Security, Audit, E-Signature) module is an optional component of the 3500 Series Data Collection Software. The SAE module provides the following functionality:

• System security – Controls user access to the software. A default Administrator user account is provided, and additional user accounts and permissions can be user-defined.

System security can be enabled or disabled globally.

- Auditing Tracks changes made to library items, actions performed by users, and changes to the SAE settings. The software automatically audits some actions silently. You can select other items for auditing and specify the audit mode. Provides reports for audited library items, SAE changes, and actions. Auditing can be enabled or disabled globally and by record type. It is enabled globally by default.
- Electronic signature (e-sig) Determines if users are permitted, prompted, or required to provide a user name and password when performing certain functions. Can be configured so that a predefined list of functions can be performed only if the data used for the functions is signed (for example, you can run a plate only if the calibration data for the system has been signed. Can be configured to require multiple signatures and to require specific users or users with specific permissions to sign.

Electronic signature can be enabled or disabled globally and by e-sig type. It is enabled globally by default.

#### Example applications

You can configure the SAE module in a variety of ways:

- Require users to log in, and leave audit and e-sig disabled.
- Allow only certain users to create or modify protocols.
- Allow only certain users to approve reviewed samples.

# Configure the security system

#### Access the Security screen and enable or disable security

The Security screen allows you to disable and enable security, control restrictions and security policies for all user accounts, and set up notifications when certain security events occur.

Security is enabled by default.

**IMPORTANT!** If you disable security, you inactivate audit and electronic signature functions. However, when you disable security, no audit record is generated to indicate that audit and electronic signature functions are disabled.

- 1. Access the Security screen.
- 2. Click **Disable** or **Enable** (Figure 34 on page 198). Note the following:
  - Disabling Security inactivates Auditing and E-signature.
  - The Disable and Enable commands are grayed when a run is in process.



- The software requires you to enter your user name and password when you enable security.
- When security is disabled, the 🕜 is not active in lower parts of the screen.

Settings Resources	🔒 Enable System Security 💣 Disable System Securi	v 😧		
AB Applied Biosystems	Account Setup			
🌮 Manage Reports			User Password	
Audit Reports	User Name		User Password	
E-Signature Reports	The length of user names must be between 0 an	d 32 characters.	The length of user passwords must be between	n 0 and 32 characters.
¥ Manage Users	Define name spacing		Define password spacing	
Users	Leading Trailing Consecutive		Leading Trailing Consecutive	
🕆 Manage Settings	Define name characteristics		Define password characteristics	
Security	Alpha Rumeric Ruppercase Lower	rcase 🗹 Special	0 Alpha 0 Numeric 0 U	ppercase 0 Lowercase 0 Special
Audi E-Signature			User may not reuse the previous 3 pass	words.
	Security Policies			6
📄 Import	Securicy Policies			
🖉 Export	Password Expiration	Account Suspensi	00	Session Timeout
Main Workflow	Passwords will expire @ tes No	Login attend	ts with an incorrect password will	Liser sessions will be timed out if
and a	every 60 days.		ts with an incorrect password will suspend the user account ○ tres	User sessions will be timed out if there is no user activity • Yes
-	Notify the user 3 days before expir		for the next 24 Hours 💌	for 60 mir
			if consecutively failing 3 time(s)	(An instrument run is not considered user activ
			within any 60 minute	

Figure 34 Security – disable or enable

## Set account setup and security policies

Security policies apply to all user accounts.

1. Under Account Setup, specify user name limits.

Account Setup	User Password
User Name	The length of user passwords must be between 8 and 32 characters.
The length of user names must be between $32$ and $32$ characters.	Define password spacing
Define name spacing	Leading V Trailing V Consecutive
Leading Trailing Consecutive	Define password characteristics
Define name characteristics	0 Alpha 0 Numeric 0 Uppercase 0 Lowercase 0 Special
Alpha Vumeric Vuppercase Vuowercase Special	User may not reuse the previous 3 passwords.

**IMPORTANT!** The software allows spaces in user names (Define name spacing). Use spaces in user names with caution. For information, see "Spaces in user names" on page 200.

- 2. Specify the *allowed* characters in user names: spaces and alpha, numeric, upper/lower case, and special characters (commas, periods, semicolons, dashes, underscores, and tildes).
- 3. Specify password limits.
- 4. Specify the *required* characters in passwords: spaces and alpha, numeric, upper/lower case, and special characters (any non-space, non-alpha, or non-numeric characters).
- 5. Specify password reuse. You cannot disable the password reuse restriction.
- 6. Under Security Policies, specify password expiration, account suspension, and session timeout settings.

Password Expiration	Account Suspension	Session Timeout
Passwords will expire  Yes No every 60 days. Notify user 3 days before expiration.	Login attempts with an incorrect password will Ves No suspend the user account Ves No for the next 24 Hours if consecutively failing 3 time(s) within any 60 minute	User sessions will be timed out if there is no user activity  Yes for 60 minutes. (An instrument run is not considered user activity.)

**Note:** A session times out while a run is in progress if the timeout period is exceeded and there is no other user activity.

- 7. Click **Setup Messaging Notification Settings** to specify when and how to notify the administrator of certain security events. For information, see "Set up messaging notifications" on page 200.
- 8. Click Save Settings.

The new settings are applied to the logged-in user the next time the user logs in.

**Spaces in user** If you allow spaces in user names, be aware of the following issues: names

- Leading and trailing spaces in user names are difficult to detect on the screen or in printed reports.
- The number of consecutive spaces in a user name is difficult to determine on the screen or in printed reports.

Spaces in user names may cause confusion when searching for an audit or E-Sig record associated with a user name. To find a record associated with a user name, you must specify the user name exactly, including leading, consecutive, and trailing spaces.

## Set up messaging notifications

1. In the Security screen (Figure 34 on page 198), click **Messaging Notifications** to display the Setup Notifications dialog box.

U	Setup Notification		
Se	t Up Notifications		
			0
	Event Name	Pop-up dialog	Message when Admin logs in
1	# of Failed Authentications over specified Time interval		
2	Session Timeout For a User		
3	Account Suspension For Failed Authentication		
4	Notification For SAE Activation		
C	Close		OK

- 2. Select the events for notification:
  - **# failed authentications over specified time interval** A user attempts to log in with an incorrect password. The message indicates the number of failed authentications.
  - Session timeout for a user No activity occurred in a user account for the specified period of inactivity.
  - Account suspension for failed authentication The user exceeds maximum number of allowed failed authentications (login attempts with an incorrect password).
  - Notification for SAE activation Security has been enabled or disabled.

- 3. Select the notification method:
  - **Pop-up dialog** The software immediately displays a pop-up message to the current user if an event is triggered by the current user. The message instructs the user to inform a system administrator of the event.
  - Message when Admin logs in If an event triggers notification, the next time any user with an Administrator role logs in, the software displays a list those events, indicating the time each event occurred and the user who triggered the event.

The Administrator has the option of acknowledging the event, which removes it from the notification list.

4. Click OK.

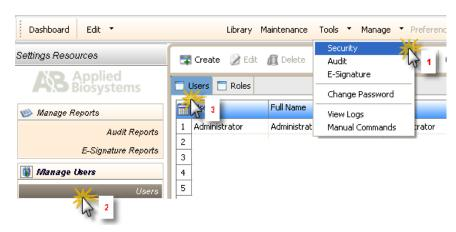
# Manage user accounts

## Create or edit a user account

The software includes a default Administrator user account with permissions (defined by the account user role) to perform all functions in the software. You cannot modify this account.

#### Create a user account

1. Access the Users screen.



2. Click **Create** to display the New User dialog box.

블 New User:					
Setup a User (2) User first name: Bi	ath Llook first appo	and Licor la	st name can not be	ometru	
	our oser hischame		st halle carrie be	empty.	
* User Name:			Created By Admin	On:	Last Updated On:
* Password:			🗹 Pre-expired		
Re-Enter Password:			Password Expires (	Dn:	
* First Name:		MI		* Last Name:	
* User Role:	Administrator 💌	* Status:	Active 🔽	Electronic Signature	🔵 Enable 💿 Disable
Email:		Phone:			
Comment	5:				< >
Close					Save

3. Enter user name, password, first name, middle initial (optional) and last name. Click a field to display the field limits, which are specified in Security settings.

**Note:** First name, MI (middle initial), and last name are used to create User Full Name, which is displayed as the name of the logged-in user.

Note: You cannot change the user name after you save the user account.

- 4. Select **Pre-expired** to require the user account to specify a new password at first log in. The Password Expires On date is specified in Security settings.
- 5. Select the user role (described in "Create or edit a user role" on page 203) and the electronic signature state (determines if a user account has permission to electronically sign objects). Leave the status set to Active.
- 6. (Optional) Enter email (for information only), phone, and comments.
- 7. Click Save.

If the Save button is grayed, it indicates an invalid entry in a field. Click a field to display the limits for the field, then enter a valid entry.

The Users screen displays the following information for each user account:

- Full Name
- Role
- Status
- Password Expired (true=yes, false=no)
- Last Modified On

- Password Change Date (by either user or administrator)
- Email (for records only)
- Phone
- Comments

# Edit a user 1. In the Users screen, select a user account, then click *B* Edit.

Note: If you select multiple users, only Status and Role will be changed.

Ē	🔄 Create 🖉 Edit 🚛 Delete 📄 View Report 🍓 Print 🝷 🕜										
	Users	Roles 6	F								
	User	Full Name	Role	Status	Password Expired	Last Modified On	Created Date	Password Change Date	Email	Phone	Comment
1	Adm	Adminis	Ad	Active	false						
2	Liser 1	User 1	Ad	Active	true	12-Feb-2009 1	12-Feb-200	12-Feb-2009 12:28:			

- 2. Edit settings as needed. You cannot edit the user name of an existing user.
- 3. Click Save.

Activate a suspended user account

2. Click **Edit**.

1. Select the user.

3. Change the status from Suspended to Active.

**Delete (inactivate)** You cannot delete a user because user records are required for auditing. To disable a user account user account, inactivate it.

- 1. Select the user.
- 2. Click **Edit**.
- 3. Change the status from active to inactive.
- 4. Click Save.

## Determine the name of the logged-in user

To display the full name of the logged-in user, place the mouse pointer on the Logout menu. The full name of the logged-in user is also displayed in the Load Plates for Run screen and the Monitor Run screen.

je	Ŧ	Preferences	Help	·	
	Æ	Administrator is	current	:ly I	ogged on.

#### Create or edit a user role

User roles determine the permissions associated with a user account.

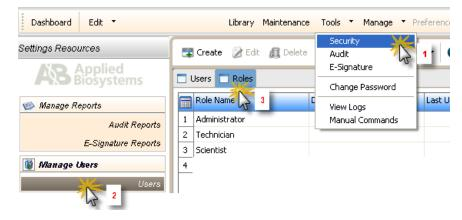
Three default user roles are included in the software. You can modify two of them, and can create your own roles with customized settings as needed:

- Administrator (cannot be edited or deleted)
- Scientist
- Technician

To determine the permissions for these roles or to edit these roles, select the role, then click  $\ge$  Edit.

Create a user role

1. Access the Roles screen.



- 2. Click **Create**.
- 3. Enter a role name and (optional) comment.
- 4. Select permissions (see Table 22 on page 204). To select all permissions in a category, select the checkbox next to the category.
- 5. Click Save Role.

#### Table 22 User role permissions

Category	Permissions
Setup	Create plate/plate template
Run	<ul> <li>Edit default instrument name</li> <li>Manage injection list</li> <li>Duplicate injection</li> <li>Re-inject</li> </ul>
Primary Analysis	Edit sample names Export sequencing results
<ul> <li>Assay</li> <li>File name convention</li> <li>Results group</li> <li>Instrument protocol</li> <li>PA protocol (primary analysis: basecalling and sizecalling)</li> <li>SA protocol (secondary analysis: sequencing, fragment analysis, HID analysis)</li> <li>QC protocol (primary analysis: HID)</li> <li>Size standard</li> <li>Dye set</li> </ul>	<ul> <li>Create</li> <li>Edit</li> <li>Delete</li> <li>Import</li> <li>Export</li> </ul>

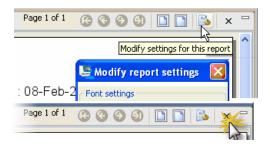
Category	Permissions
Plates and templates	<ul> <li>Edit</li> <li>Delete</li> <li>Import</li> <li>Export</li> </ul>
Locking/Unlocking	<ul> <li>Assay</li> <li>File name convention</li> <li>Results group</li> <li>Instrument protocol</li> <li>PA protocol</li> <li>SA protocol</li> <li>QC protocol</li> <li>Size standard</li> <li>Dye set</li> </ul>
Preferences	<ul> <li>Edit system preferences</li> <li>Export system preferences</li> <li>Import system preferences</li> <li>Export user preferences (all)</li> </ul>
Calibrations	<ul><li>Perform spatial calibration</li><li>Perform spectral calibration</li></ul>
Performance check	Run performance check install standards
Archiving	<ul><li>Archive</li><li>Purge</li><li>Restore</li></ul>
SAE configuration	<ul><li>Configure SAE</li><li>Log in to timed-out user sessions</li></ul>

Table 22 User role permissions (continued)

- Edit a user role 1. In the Roles screen, select a user role, then click *P* Edit.
  - 2. Edit settings as needed. You cannot edit the Administrator user role.
  - 3. Click Save.

#### View and print a user report

- 1. Select the User or Roles tab. Click 📄 View Report.
- 2. In the Report screen, click toolbar options to manipulate the report as needed. Place the mouse pointer over an item for a description of the item.
- To print the report, click Print. Close the report.



Save electronic To copies (.pdf) of the report

To save the report electronically (.pdf), print the report and select **CutePDF Writer** as the printer.

## Example reports

User	Report				g						
#	User	Full Name	Role	Email	Phone	Status	Created Date	Last Modified On	Password Change Date	Password Expired	Comments
1	Administra tor	a Administr ator	Adminis trator			A ctive				false	
2	technician	First Name MI User2	Technici an			A ctive	29-Jan- 2009 10: 13:57 A M	01-Feb- 2009 10: 10:49 AM	31-Jan- 2009 10:12: 48 A M	false	
3	scientist	FirstNam e MI User3	Scientist			A ctive	29-Jan- 2009 11: 30:12 A M	01-Feb- 2009 10: 10:22 AM	31-Jan- 2009 03:36: 58 P M	false	
4	Analyst1	First Name MI Analyst1	Log in to timed- out scssion			A ctive	01-Feb- 2009 10: 21:26 A M	01-Feb- 2009 10: 21:26 AM	01-Feb- 2009 10:21: 26 A M	true	
	User R	ole Rep	ort								
	#	Rol	e Name		De	scriptio	n	Last Up	dated Date	Use	r Counts
	1	Adm	inistrator								3
	2	Tec	hnician								1
	3	Sc	ientist								1
	4 L	_og in to tin	ned-out se	ssion				01-Feb-200	9 10:12:18 AN	1	1

Tools 🔻 Manage 🝸 Preference:

Security

E-Signature Change Password

View Logs

Manual Commands

Audit

# Manage auditing

## Access the Audit screen and enable or disable auditing

The Audit screen controls the auditing state (enabled/disabled), the events that are audited, and the reasons available to users when audit mode is set to Prompt or Required.

Auditing is enabled by default.

**IMPORTANT!** If you disable security, you inactivate audit and electronic signature functions. No audit record is generated for the inactivation of audit and electronic signature functions when you disable security.

Maintenance

- 1. Access the Audit screen.
- 2. Click **Disable** or **Enable** (Figure 35 on page 207).

**Note:** When auditing is disabled, the **?** is not active in lower parts of the screen.

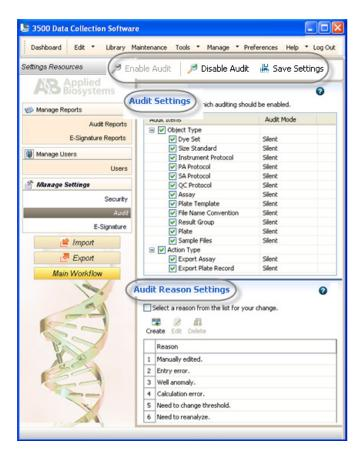


Figure 35 Audit – disable or enable

# Select objects to audit

Object Type (audit records displayed in Object Audit History)	Action Type (audit records displayed in Action Log)
Dye set	Export assay
Size standard	Export plate record
Instrument protocol	
<ul> <li>PA protocol (primary analysis)</li> </ul>	
SA protocol (secondary analysis)	
QC protocol	
Assay	
Plate template	
File name convention	
Results group	
Plate	
Sample files	

1. Select the objects and actions to audit and the mode for each enabled item.

**Note:** For a list of items that the system audits silently in addition to the configurable items listed above, see "Generate audit reports" on page 209.

- 2. Set the Audit Mode for each item you enable for auditing:
  - **Prompt** 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.
  - Silent The event is audited, no reason prompt is displayed.
- 3. Click Save Settings.

## Create audit reason settings

You can create, modify and delete the reasons that are available for selection in the Audit Reason dialog box (displayed when a user performs an audited action).

1. To require users to Audit Reason Settings select a pre-defined reason in the Audit Select a reason from the list for your change. Reason dialog box 廁 Create Edit Delete 🔚 Audit Reason (displayed when a user performs an Setup Audit Reason Reason audited action), 🐼 Audit Reason is a required field. Provide a un 1 Manually edited. value. enable the User must 2 Entry error. 3 Well anomaly. select a reason Enter an Audit Reason in t 4 Calculation error. checkbox. Users are 5 Need to change threshold. not permitted to enter 6 Need to reanalyze. a reason. Close 2. As needed, click 🔜

Create, or select a reason, then click 📑 Edit or 慮 Delete.

# Generate audit reports

# **Display audit histories**

1. Access the Audit Reports screen.

**Note:** To access the Audit Reports screen, the user role for an account must specify the Configure SAE permission. Users without the Configure SAE permission can view object audit histories for individual entries in the libraries by selecting entries, then clicking View Audit History (see "View audit and e-signature histories for library entries" on page 142.



- 2. Select a tab to display:
  - **Object Audit History** The most recent audit for all user objects (samples and objects in the Library) that have been audited.
  - System Configuration History SAE configuration records, including audit history for each user account.
  - Action log System-specified audit events.
- 3. (Optional):
  - Sort the table. See "Multi-column sorting" on page 72.
  - Specify filters (date range, user name, action, object or record type, object or record name, reason), then click **Go**.

Note: The Reason field in System Configuration History is not used.

- Select a record, then click **Show Object History** or **Show Audit Details**.
- In the history dialog box, select a record, then click **Show Audit Details**.
- Click **Table Settings**, then specify the columns to show or hide.

# Review the object audit history

Audit records The Object Audit History lists the most recent audit for the user objects listed below (samples and objects in the libraries) that have been audited.

- Dye set
- Size standard
- Instrument protocol
- PA protocol (primary analysis)
- SA protocol (secondary analysis)
- QC protocol

- Assay
- Plate template
- File name convention
- Results group
- Plate
- Sample files
- Audit action Possible actions for all objects are update, create, and delete. Audit records are generated under the following conditions:

Action	Description
Update	The auditing of updates depends on whether an object is modified or overwritten:
	• Modified – A record is created when an object is modified.
	• <b>Updated</b> – A record is not created when an object is overwritten in the library. Example: You create a plate, then create a results group from within the plate and save it to the library. You then open the plate, edit the results group from within the plate, then save it to the library. A message indicates that the results group already exists and asks if you want to overwrite it. You click Yes. This action is considered a creation of a new results group, not a modification of the existing results group No Update record is created; a Create record is created.
Create	A record is created when you:
	Create an item in the library.
	Create an item from within another item.
	<ul> <li>Modify an item from within another item, then overwrite the item in the library when you save it (as described in the "Updated" bullet above).</li> </ul>
Delete	The auditing of deletions depends on the item deleted:
	• <b>Items in the library</b> – A record is retained until it is deleted from the library. The deletion of the item from the library is not audited. For example, if you <i>delete</i> a size standard from the library, no audit record for the deletion is listed in the Object Audit History.
	• Items within other items – The deletion of an item from within another item is audited. For example, if you <i>change</i> the size standard in a QC protocol, an audit record for the change (considered a deletion) is listed in the Object Audit History.



# Display the<br/>object historyTo display the history for an object, select the object, then click Show Object<br/>History.

The object history shows the audit history for the object and for all objects contained in the selected object. For example, when you create an assay, a copy of the instrument protocol, the primary analysis protocol (and therefore dye set, and size standard), and the secondary analysis protocol are included in the assay object. The objects contained within an object have audit histories distinct from the audit history of the objects stored in the Library.

# Review the system configuration history

The System Configuration History lists SAE configuration records.

Note: The Reason field in System Configuration History is not used.

Record Type	Action	Corresponds To
Security settings	Update	<ul> <li>Enable security</li> <li>Disable security</li> <li>Modify security policies:</li> <li>Session timeout settings</li> </ul>
Account settings	Update	<ul> <li>Modify user name settings</li> <li>Modify password settings</li> <li>Modify security policies:</li> <li>Password expiration</li> <li>Account suspension</li> </ul>
Audit reason for change	Update	Modify reason for change
	Create	Create reason for change
	Delete	Delete reason for change
Audit settings	Update	Enable auditing Disable auditing
Audit type	Update	Modify audit settings
Audit type	Update	<ul><li>Modify audit settings</li><li>Create reasons for change</li><li>Delete reasons for change</li></ul>
E-Signature function	Update	<ul> <li>Modify the number of signatures or the authorities for a "prompt before" function</li> <li>Modify the Enable state of either a "check before" or "prompt before" function</li> </ul>
E-Signature settings	Update	<ul><li>Enable e-signature</li><li>Disable e-signature</li></ul>
E-Signature type	Update	<ul><li>Modify e-signature settings</li><li>Modify the enable state of an E-Signature Type</li></ul>

#### Table 23 Audit – system configuration history

Record Type (continued)	Action	Corresponds To
Role assignment	Create	<ul> <li>Create a new user account</li> <li>Assign a different user role to an existing user account</li> </ul>
	Delete	Assign a different user role to an existing user account
Role permissions	Update	Modify user role permissions
	Create	Create a user role - creates one role assignment record for each permission in a role
	Delete	Delete a user role - creates one role delete record for each permission in the deleted role
User account	Update	<ul><li>Edit</li><li>Suspend</li></ul>
	Create	Create new user account
User role	Update	Modify user role
	Create	Create user role
	Delete	Delete user role

Table 23	Audit – system	configuration	history	(continued)
----------	----------------	---------------	---------	-------------

# Review the action log

The Action log lists system-specified audit events.

All items in the action log are audited silently, except for the items noted as configurable. Configurable items may include comments in the action log.

Table 24	Audit – action log
----------	--------------------

Category	Action
Assay	Assay exported successfully
	<b>Note:</b> Only one audit record is generated if you export multiple assays.
Log In	<ul><li>User logged in</li><li>Login failed</li><li>User logged out</li></ul>
Maintenance Wizards	<ul> <li>Remove Bubbles Wizard started</li> <li>Flush Array Port Wizard started</li> <li>Change Polymer Type Wizard started</li> <li>Change Array Wizard started</li> <li>Replenish Polymer Wizard started</li> <li>Perform Fill Polymer Wizard</li> <li>Perform Water Wash Wizard</li> </ul>
Plate	Plate exported successfully <b>Note:</b> Only one audit record is generated if you export multiple plates.

Category	Action
Run	<ul> <li>Start</li> <li>Pause</li> <li>Resume</li> <li>Stop (Abort injection)</li> <li>Terminate (injection list)</li> </ul>
SAE Configuration	Export
System Audit Records	<ul><li>Archive</li><li>Purge</li><li>Restore</li></ul>
System Action Records	<ul><li>Archive</li><li>Purge</li><li>Restore</li></ul>
User Profile	Export

Table 24 Audit – action lo	oq (continued)
----------------------------	----------------

## View and print audit reports

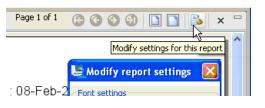
- 1. Display the records of interest as described above.
- 2. Filter the list to decrease the time required to generate reports.

**IMPORTANT!** You cannot cancel a report after you click a view button.

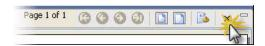
3. Click View Audit Summary Report or View Audit Detailed Report.

Sy	/ste	m C	onfig	juration	History Sum	nm ary	/ Report				
	#		Date		User Name	User	Full Name	Recor	d Type	Record Name	Action
	1	28-Jar	n-2009 ( PM	05:01:08	A dministrator	Ad	Administrator Secu		ecurity Settings		Update
	2	28-Jar	n-2009 ( PM	05:00:57	A dministrator	Ad	ministrator	Security	Settings		Update
Sys	ster	n Co	nfigu	uration H	istory Detai	led R	eport				
1	Dat	e:		28-Jan-200	9 05:01:08 PM		Action :	U	pdate		
	User Name : Administrator			User Full Nar	ne: A	dministra	tor				
	Rec	ord Ty	ype:	Security Se	tings		Record Nam	e:			
	Ħ	ŧ	Reco	rd Type	Object Na	me	Old Va	lue	Cu	rrent Value	Action
	1		Securi	ty Settings	Security On / Se Off	ecurity	DISABI	ED	E	ENABLED	Update
2	Dat	e:		28-Jan-200	9 05:00:57 PM		Action :	U	Ipdate		
	Use	r Nam	ie :	Administrat	or		User Full Nar	ne: A	dministra	tor	
	Rec	ord Ty	ype:	Security Se	ttings		Record Nam	e:			
	Ħ	ŧ	Reco	rd Type	Object Na	me	Old Va	lue	Cu	rrent Value	Action
	1		Securi	ty Settings	Security On / Se Off	ecurity	ENABL	ED	C	DISABLED	Update

4. In the Report screen, click toolbar options to manipulate the report as needed. Place the mouse pointer over an item for a description of the item.



- 5. To print the report, click **Print**.
- 6. To save the report electronically (.pdf), print the report and select **CutePDF Writer** as the printer.
- 7. Close the report.



#### Archive, purge, and restore audit records

The audit archive function makes a copy of audit records. Purge makes a copy of audit records, and then deletes them. You can use the Restore function to restore purged audit records.

For information on archiving library items (datastore), see "Archive, purge, and restore data" on page 254.

Archive and	To selectively archive or purge (delete) system configuration or action audit records:
purae	

1. Select records in the appropriate screen.

- 2. Click 📩 Archive Audit Records or 🕮 Purge Audit Records.
- 3. If you select Archive, specify a location and name for the .asz archive file.
- **Restore** To restore system configuration or action audit records, click **Restore**, then select the .asz file to restore.

## Export audit records

As needed, you can export audit records to a .txt file for additional manipulation and reporting outside the 3500 Series Data Collection Software.

- 1. Display the records of interest as described above.
- 2. Click 🛃 Export Audit Records.
- 3. Specify a name and location for the export .txt file.
- 4. Click Save.

**Note:** If you export audit records for samples that are not in their original location (samples have been deleted or moved), an error message is displayed. Return sample data files to their original location, then export again.

# Manage electronic signature

## Access the E-Signature Settings screen and enable or disable e-sig

**IMPORTANT!** If you disable security, you inactivate audit and electronic signature functions. No audit record is generated for the disabling of audit and electronic signature functions when you disable security.

- 1. Access the E-Signature Settings screen screen.
- 2. Click **Disable** or **Enable** (Figure 36 on page 216).

**Note:** When e-sig is disabled, the **?** is not active in lower parts of the screen.



Dashboard Edit *	Library Maintenance	Tools * Manage * Preferen	ces Help * Log Out
Annied CA	Enable E-Sig i Disable E-Sig i B Sectors	Save Settings	
Manage Reports	E-Signature Type Sele	ect the functions after which	the system will pror
Audit Reports E-Signature Reports	Approve Dye Set Approve Size Standard Approve Spatial Calibration	Function	
Manage Users	Approve Spectral Calibration     Approve Instrument Protocol		
Users	Approve Instrument Protocol Approve Sizecall Protocol Approve Basecall Protocol		
Manage Settings	Approve QC Protocol		
Security Audit	Approve Gene Mapper Protocol     Approve Gene Mapper IDx Protocol     Approve SeqScape Protocol     Approve MicroSeg ID Protocol		
E-Signature	Approve Assay	4	2
ピ Import	L Hyprovernake	ect the functions before whi	ch the system will ch
Export	Approve Sample Approve Sequencing Install Standard Results	Function	# Signatur
Main Workflow	Approve MicroSeqID Install Standard Results     Approve Fragment Install Standard Results     Approve HID Install Standard Results		
10		S	

Figure 36 E-Sig – disable or enable

## Select the actions that allow signature

**IMPORTANT!** Do not change electronic signature settings during a spectral calibration.

1. Select the E-Signature Settings checkbox next to Select the electronic signature types that should be allowed. an item in the Select the functions after which E-Signature Type **E-Signature** Type Approve Dye Se Function Approve Size Standard list to identify Save Dve Set Approve Spatial Calibration events for which Approve Spectral Calibration to allow

electronic signature (see Table 25 on page 218). This selection activates the E-Sig button for the selected items; it does require an electronic signature for these selections.

- 2. (Optional) For each item that you select,:
  - a. From the top-right of the screen, select a function *after which* the system will prompt for electronic signature. This selection presents an e-sig prompt to users when they perform a function. Users can sign or can continue without signing.
  - b. From the bottom-right of the screen, select a function (start run) before which the system will check for required electronic signatures (see Table 26 on page 219). This selection presents an e-sig prompt to users when they start a run if the required signatures have not previously been made. Users must sign before they can continue. For "check before" functions, you can also:
    - Change the number of signatures required.
    - Set a special authority for a signature: click the Authorities Required field, then select the user account or the user role to require for electronic signature of this function. By default, each required signature needs no special authority; any user can sign.
    - Click Apply.

	E-Signature Type		Select the rund	tions after which the systi	em will prompt for an electro	nic signatui
ł	🖌 🗹 Approve Dye Se	et	Function			
L	Approve Size St	andard		and an Information		
	Approve Spatia	l Calibration 🛛 💦 🦳	Accept Sp	atial Calibration		
l	Approve Spectr	al Calibration				
l	Approve Instru	ment Protocol	Select the func	tions before which the sys	stem will check for required (	electropic si
	Approve Sizeca	ll Protocol	boloce the rane	dons berore which are sy.	com will check for required (	Sideer of the St
	Approve Baseca	all Protocol	Function	# Signatures Required	Authorities Required	
	Approve QC Pro	otocol	Start Run	-	· · · · · · · · · · · · · · · · · · ·	
	🗹 Approve Gene I	Mapper Protocol			Any	
l	🗹 Approve Gene I	Mapper IDx Protocol	1			
	Approve SeqSc	ape Protocol	NE			
l	Approve MicroS	ieq ID Protocol	-Select the mi	nimum signatures that mus	t exist for data to qualify a	s heina sian
	🗹 Approve Assay		Delect the fill	ninan signacaros criacina.	a conserver adea co quality a	s boing sign
	🗹 Approve Plate 1	ſemplate	Function Na	me: Run.Start		
	🗹 Approve Plate					
	🗹 Approve Sample	э	# Signature	Doguirod	1	
	Approve Seque	ncing Install Standard Results	* Signature:			
	Approve MicroS	eqID Install Standard Results				
		ent Install Standard Results	Authorities	Required Any		
	Approve Fragm					
		stall Standard Results				
		stall Standard Results				

3. Click Save Settings.

By default, no E-Signature types are enabled.

Table 25	E-signature settings to prompt after
----------	--------------------------------------

E-Signature Type	Function to Prompt After
Approve Dye Set	Save
Approve Size Standard	Save
Approve Spatial Calibration	Accept
Approve Spectral Calibration	Accept
Approve Instrument Protocol	Save
Approve Sizecall Protocol	Save
Approve Basecall Protocol	Save
Approve QC Protocol	Save
Approve Size Standard	Save
Approve Spatial Calibration	Accept
Approve Spectral Calibration	Accept
Approve Instrument Protocol	Save
Approve Sizecall Protocol	Save
Approve Basecall Protocol	Save
Approve QC Protocol	Save
Approve GeneMapper Protocol	Save
Approve GeneMapper IDX Protocol	Save
Approve SeqScape Protocol	Save

E-Signature Type	Function to Prompt After
Approve MicroSeq ID Protocol	Save
Approve Assay	Save
Approve Plate Template	Save
Approve Plate	Save
Approve Sample	Save
Approve Sequencing Install Standard Results	Accept
Approve MicroSeq ID Install Standard Results	Accept
Approve Fragment Install Standard Results	Accept
Approve HID Install Standard Results	Accept

Table 25 E-signature settings to prompt after (continued)

#### Table 26 E-signature settings to check before

E-Signature Type	Function to Check Before	Signatures and Authorities Required (defaults if enabled)
Approve Spatial Calibration	Start Run	1 signature, any authorities
Approve Spectral Calibration	-	(any user, any user role)
Approve Spatial Calibration	-	
Approve Spectral Calibration	-	
Approve Plate	-	
Approve Sequencing Install Standard Results	_	
Approve MicroSeq ID Install Standard Results	_	
Approve Fragment Install Standard Results	-	
Approve HID Install Standard Results	-	

#### How the software prompts electronic signature before a run

If the system is configured to check that data is signed fore starting a run and the data for the run is not signed, a message is displayed when the user clicks **Start Run**.

#### Example

The e-sig system is configured to require signatures from two users (one from the user account named Administrator, and the other from any user account with a scientist user role) for a spatial calibration before it can be used in a run. The spatial calibration has not been signed.

A user starts a run. The following message is displayed:

🕒 E-Si	gnature Check
?	The Spatial Calibration Spatial_Run 2009-01-29-09-59-13 record must be signed by 2 users to be utilized for this Start Run, but is not signed by any users. The following signatures are missing: • Approve Spatial Calibration by a user with User Account: Administrator authority • Approve Spatial Calibration by a user with User Role: Scientist authority Would you like to sign now?
	Yes No

Before the run can start, the following users must sign:

- The Administrator user
- Any other user with the Scientist role specified and electronic signature enabled in their user account

If a user that does not meet the specified criteria signs, this message is displayed again.

# Generate e-signature reports

# Display e-signature records

- 1. Access the E-Signature Reports screen.
- 2. (Optional):
  - Specify filters (date range, user name, action, object type, object name), then click **Go**.
  - Select a record, then click Show Object History.
  - In the history dialog box, select a record, then click **Show E-Signature Details**.
  - Double-click column headers to sort. Multi-column sorting is supported (see "Multi-column sorting" on page 72).



- Customize the table (see "Customize tables" on page 72).
- 3. The records that are displayed (if they are specified in E-Signature settings) are:
  - Approve Dye Set
  - Approve Size Standard
  - Approve Spatial Calibration
  - Approve Spectral Calibration
  - Approve Instrument Protocol
  - Approve Sizecall Protocol
  - Approve Basecall Protocol
  - Approve Qc Protocol
  - Approve Genemapper Protocol
  - Approve Genemapper ID-X Protocol
  - Approve Seqscape Protocol

- Approve Microseq ID Protocol
- Approve Assay
- Approve Plate Template
- Approve Plate
- Approve Sample
- Approve Sequencing Install Standard Results
- Approve Microseq ID Install Standard Results
- Approve HID Install Standard Results
- Approve HID Install Standard Results

## View and print e-signature reports

1. Display the records of interest as described above.

Note: Filter the list to decrease the time required to generate reports.

2. Click View E-Sig Summary Report or View E-Sig Detailed Report.

E-Signature Summary Report					lay-2009 03:13:09 PM Software Version 1.0.0	
#	Date	User Name	User Full Name	Object Type	Object Name	Comments
1	04-May-2009 03:11: 47 PM	Administrator	Administrator	Approve Spatial Calibration	Spatial_Run 2009-05-04-15- 10-51	Spatial calibration is acceptable



E-Signature Detailed Report

User Name : Administrator

User Name : Administrator	User Full Name : Administrator
Object Type : Approve Spatial Calibration	Object Name : Spatial_Run 2009-05-04-15-10-51
Date : 04-May-2009 03:11:47 PM	Comments : Spatial calibration is acceptable
Esignature Details Esignature Type : Approve Spatial Calibration	
Signed By : Administrator	
Full Name : Administrator	
Signed On :04-May-2009 03:11:47 PM	
Authority : User Account: Administrator,	User Role: Administrator

Object Details

5	Intensity	16 5 -5 5 13 2 8 9 12 -1 -1 7 3 6 -1 1 -7 4 4 8 5 -2 0 9 2 -3 0 -6 -6 -2 7 -4 -4 -9 -3 -8 -3 8 -7 4 2 1 -3 -5 -2 1 -3 -4 5 6 5 3 0 2 0 1 -1 2 5 1 3 2 3 0 1 3 0 1 -3 -4 0 -5 -7 -5 -7 -3 1 4
6	Run ID	Spatial_Run 2009-05-04-15-10-51
7	Number of Capillaries	24
8	Locked	false
9	Instrument	13527-029
10	Capillary Array	80K0850

- 3. In the Report screen, click toolbar options to manipulate the report as needed. Place the mouse pointer over an item for a description of the item.
- 4. To print the report, click **Print**.
- 5. To save the report electronically (.pdf), print the report and select CutePDF Writer as the printer.

Page 1 of 1	6 6 9 9 🗅 🗋 🔖 ×	
	Modify settings for this report	^
	🖶 Modify report settings 🛛 🔀	
:08-Feb-2	Font settings	
	Select the font to be used in reports.	
	10 🗸	

Report created on : 04-May-2009 03:14:16 PM 3500 Data Collection Software Version 1.0.0 6. Close the report.



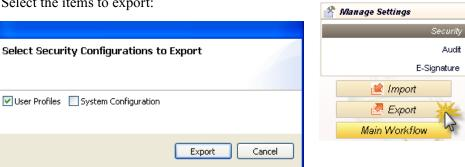
# Export e-sig records

As needed, you can export e-sig records to a .txt file for additional manipulation and reporting outside the 3500 Series Data Collection Software.

- 1. Display the records of interest as described above.
- 2. Click **Export E-Sig Records**.
- 3. Specify a name and location for the export .txt file.
- 4. Click Save.

# Export and import user accounts, security, audit, and electronic signature settings

- 1. In any screen in the SAE module, click 🛃 Export Export in the navigation pane.
  - 2. Select the items to export:



- User Profiles Contains all settings in the following screens:
  - Edit User All user accounts with Active status
  - User Role All user roles and associated permissions (in case a user account specifies a user role that does not exist on the system into which you import the profiles)
- System Configuration Contains all settings in the following screens:
  - Security Account setup and security policies
  - Audit Objects selected for auditing, audit modes, and reasons
  - E-Signature Settings Objects selected for E-Signature, functions, number of signatures, and authorities
  - User Roles All user roles and associated permissions
- 3. Click Export.
- 4. Specify the name and location for the exported .dat file, then click Save. A message is displayed when the export completes.
- Import 1. In any screen in the SAE module, click **Import** in the navigation pane.
  - 2. Select the .dat file to import, then click **Open**. A message is displayed asking if you want to overwrite the current system configuration. Click Yes.

If any imported user accounts already exist on the system, you are prompted to overwrite or skip each account.

Please resolve conflicts				
	User	Overwrite	Skip	
1	Administrator			
2	User 1			

🧃 Manage Users

Users

# Section 2 Users

# Users overview of System Security, Audit Trail and E-Signature

The Security, Audit, E-Signature (SAE) module is an optional component of the 3500 Series Data Collection Software. The SAE module provides the following functionality:

- System security Controls user access to the software.
- Auditing Tracks changes made to library items, actions performed by users, and changes to the SAE settings.
- Electronic signature (e-sig) Requires users to provide a user name and password when performing certain functions.

Depending on the way that your administrator configures these features, you may see the following dialog boxes and prompts when you use the software.

# Security

Log in If security is enabled on your system, you must provide a user name and password to access the software.

Your access to functions in the software is based on the permissions associated with your user account. Functions for which you do not have permissions are grayed out.

If your system is configured for

블 3500 Log In	
3500 Log In	
Provide your user name and password to login.	
User Name:    Password:	
Reset	

password expiration, you will periodically be prompted to change your password. If your system is configured to monitor failed log in attempts, you will be locked out of the software if you incorrectly enter your user name or password for a specified number of times.

**Permissions** If your user account does not have permission to perform any function in the software, menu commands are grayed.

To display the full name of the logged-in user, place the mouse pointer on the Logout menu. The full name of the logged-in user is also displayed in the Load Plates for Run screen and the Monitor Run screen.		
When your password is about to expire, a message is displayed when you log in.		
To change your password, select <b>Tools ► Change</b> Password.  Maintenance Tools ▼ Manage ▼ Preferences Security		
Enter your current password, then enter the new password two times, then click <b>OK</b> .  Audit E-Signature  Change Password View Logs Manual Commands		
If your system is configured to suspend a user account for failed logins, and you enter an incorrect user name and password for more than the allowed number of times, your user account is suspended, and the Log In dialog box indicates that your account is inactive.		
<ul> <li>There are two ways to activate a suspended account:</li> <li>You can wait until the suspension period ends.</li> <li>An administrator can change the account status from Suspended to Active.</li> </ul> Note: While a user is suspended, another user can click <b>Reset</b> , then log in and		
replace the suspended user.		
If your system is configured to timeout and there is no user activity for the specified time, the Log In dialog box indicates that your user session has timed out. You must enter your user name and password to access the software. Note: The administrator or another user with permission to log in to timed-out sessions can click <b>Reset</b> , then log in.		

# Audit

If your system is configured for auditing, you may be prompted to specify a reason when you make certain changes in the software.

Based on your system configuration, you can either select a reason or enter a reason for change.

🔚 Audit F	leason		X
Setup Au	dit Reason		
Select from	List of Reasons:		~
Comments:		Manually edited. Entry error. Well anomaly. Calculation error.	
Close		Save	

# **Electronic signature**

If your system is configured for electronic signature, you may be prompted to provide your user name and password when you perform certain actions in the software.

If an item is set to require two signatures, the signers are not required to sign at the same time. When the first signer signs, the E-Sig status is set to Partially Signed. When the second signer signs, the E-Sig status is set to Signed.

블 Electronic Si	gnature	×
Electronic Signature Enter your user name, password and any comments.		
E-Sig Record Type:	Approve Spatial Calibratic 🐱	
User Name:	User1	
Password:		
Comments:		
	OK Cancel	

You may also be permitted to sign objects such as plates, calibrations, or other library items. If electronic signature is enabled for items, any of the following may apply:

- The 🗯 E-Signature button is enabled in the library or the calibration.
- You are prompted to sign as described in "How the software prompts electronic signature before a run" on page 220.

• The Open Plates dialog box or the library displays an "Is signed" column that reflects the electronic signature status of an item.

Open Plate From Library								
	ructions ct row from table and click	on "Open" buttor	n.					
Filter: HID Search: All So Clear @								
	Plate Name	Туре	1	te	Date Modified	(	Is Signed	
	Plate 01	HID			12-Apr-2009 05:13:13 PM	Г	No	
1								

# Maintenance schedule

This section lists the common tasks required to maintain your Applied Biosystems 3500/3500xL Genetic Analyzers in good working condition.

The Dashboard, in conjunction with the data entered in the schedule section of the Planned Maintenance, provide a comprehensive outline of maintenance tasks.

**WARNING!** Wear appropriate protection, including gloves, laboratory goggles, and coat whenever you work with the fluids used on this instrument, or parts that may come into contact with these fluids.

**IMPORTANT!** Use the cleaning agents as described in this manual, only. Use of cleaning agents not described in this manual can impair the instrument.

For the instrument troubleshooting issues, see Appendix E, "Troubleshoot" on page 299.

## **Review maintenance notifications**

Review maintenance notifications list in the Dashboard daily, then perform the scheduled tasks.

Name	Priority	Notification Date	Description	Action	
Perform Performance Check	HIGH	28-Jan-2009 12:00:00 AM	Performance Check	1	×
Clean Drip Tray	HIGH	28-Jan-2009 12:00:00 AM	Clean Drip Tray	×	x
Clean Autosampler	HIGH	28-Jan-2009 12:00:00 AM	Clean Autosampler	×	x
Replace Reservoir Septa	HIGH	28-Jan-2009 12:00:00 AM	Replace Reservoir Septa	×	x
Wash Pump Trap	HIGH	28-Jan-2009 12:00:00 AM	Wash Pump Trap	×	x

When you complete a task, click  $\bigvee$  to mark it as complete, click  $\bigotimes$  to mark it as dismissed.

**Note:** Completed and dismissed tasks are removed from the Maintenance Notification section, and they do not appear again unless they are repeating tasks. Dismissed tasks can be logged in the Notifications Log.

All actions are recorded in the Notification Log. See "Review the Maintenance Notifications Log" on page 257.

# Daily instrument maintenance tasks

Clean the assemblies, anode buffer container, and cathode buffer container, and ensure that the outside of the assemblies is dry.

**IMPORTANT!** Use the cleaning agents as described in this manual, only. Use of cleaning agents not described in this manual can impair the instrument.

Task	Frequency	For information, see
Check consumables on the Dashboard – Refer to the gauges on the Dashboard to see the status for anode buffer container, cathode buffer container, and polymer.	Before each run	"Check consumables on the Dashboard" on page 236
Visually inspect the level of fluid inside the anode buffer container and the cathode buffer container. The fluid must line up with the fill line.		"Change the cathode buffer container (CBC)" on page 238
Ensure that the plate assemblies are properly assembled.	-	"Prepare the plate
<b>IMPORTANT!</b> Align the holes in the plate retainer with the holes in the septa to avoid damaging capillary tips.		assembly" on page 53
Ensure that the plate assemblies and the cathode buffer container are positioned on the plate deck properly. They should sit securely on the deck.	-	"Load the plate in the instrument" on page 53
Ensure the array locking lever on the capillary array is secured.	-	Chapter 1, Instrument and Software Description
Check for bubbles in the pump block and channels.	Daily or before each run	"Remove bubbles from
Note: Use the Remove Bubble wizard to remove bubbles.		the polymer pump" on page 251
Check the loading-end header to ensure that the capillary tips are not crushed or damaged.	-	"To change the capillary array" on page 252
Ensure that the pump block is in pushed back position.	Daily	Chapter 1, Instrument and Software Description
Clean the instrument surfaces of dried residue, spilled buffer, or dirt.	-	"Routine instrument cleaning" on page 242
Check for leaks and dried residue around the Buffer-Pin Valve, check valve, and array locking lever.	-	"Check maintenance notifications" on
IMPORTANT! If leaks persist, contact Applied Biosystems.		page 28

# Weekly instrument maintenance tasks

Task	Frequency	For information, see
Check the storage conditions of the used arrays to ensure the array tip is covered in the reservoir.	Weekly	"Check stored capillary arrays" on page 240
Run the Wash Pump and Channels wizard.		"Wash the pump chamber and channels" on page 249
Use a lab wipe to clean the anode buffer container valve pin assembly on the polymer delivery pump.		Chapter 1, Instrument and Software Description
Restart the computer and instrument.		"Reset the instrument" on page 314.

# Monthly instrument maintenance tasks

Task	Frequency	For information, see
Flush the pump trap	Monthly or as needed	"Flush the water trap (pump trap)" on page 241
Empty the condensation container and the water trap waste container. The waste container is to the right of the pump block.		Chapter 1, Instrument and Software Description
Replace cathode buffer container septa.		"Change the cathode buffer container (CBC)" on page 238
Run a performance check	_	Chapter 5, Calibrate and Check Performance
Clean the autosampler	-	"Routine instrument
Clean the drip tray	-	cleaning" on page 242
Check disk space	-	"Monitor disk space" on page 256
Defragment the hard drive	Monthly Before fragmentation reaches 10%.	"Defragment the computer hard drive" on page 257

# Quarterly maintenance tasks

Task	Frequency	For information, see
Run performance check	Every three months	Chapter 5, Calibrate and Check Performance

## Annual planned maintenance tasks

Call your Applied Biosystems representative to schedule annual planned maintenance.

## As-needed instrument maintenance tasks

Task	Frequency	For information, see	
Change the tray.	As needed	"Routine instrument cleaning" on page 242	
Remove dried polymer from the capillary tips with a lint-free wipe moistened with deionized water.			
Archive and purge library objects		Chapter 6, Manage	
Dashboard  Manage  Archive or Purge		Library Resources	

# Use the maintenance calendar

The Maintenance calendar is a monthly or daily view of the routine maintenance tasks scheduled for your instrument. When a task is due to be performed, it is listed in the Maintenance Notifications list in the Dashboard (see "Review maintenance notifications" on page 229).

## View the calendar

To go to the Schedule from the Dashboard:

1. In the Dashboard, click Maintain Instrument toggle key.

The Planned Maintenance options appear on the left-hand pane, highlighted below:

The Dashboard provides you with a list of current maintenance notifications, as shown. Click ? for information.

- 2. From the Left-hand pane, under Planned Maintenance, click **Schedule**
- 3. Click ? on the top left-hand corner of the Schedule for more information.

Additionally, Applied Biosystems suggests that you add the regular maintenance tasks listed below to the maintenance calender.



#### **Default calendar entries**

A set of Applied Biosystems-recommended tasks are scheduled in the calendar, flagged with FR (Factory Repeating) in the monthly view and F (Factory) in the daily view. User-specified repeating tasks are flagged with R (Repeating) in the monthly view, see picture below.

You can change the priority of factory tasks, but you cannot remove them from the calendar or alter the frequency at which the notifications for the tasks are displayed.

Additionally, Applied Biosystems suggests that you add to the maintenance calendar:

- The regular maintenance tasks.
- A maintenance task to replace a consumable based on its installation date (for example, create a task to replace the polymer for two days before the polymer will expire)

#### Create calendar entries

To create a new scheduled task, click Create and follow the prompts.

The following is an example of scheduled events in the calender.



The Month and Day tabs allow you to view your schedule in different formats. Click **Detach** to move the calendar window.

Dashboard Edit •		
💥 Maintenance	🗔 Create 🛛 🖉 Edit	🕼 Delete 🛛 Detach 🛛 🚱
AB Applied Biosystems	🚱 Month 🔮 Day	
Spatial	Monday	Tuesday

# **Review the Maintenance Notifications Log**

The Notifications Log is a history of all notifications messages and the action taken for the task (completed or dismissed). You can use this option to review a previous run information.

The Dashboard provides you with a list of current routine and maintenance notifications, as explained below.

Multi-column sorting is supported (see "Multi-column sorting" on page 72.

To go to the Notifications Log from the Dashboard:

#### 1. Click Maintain Instrument

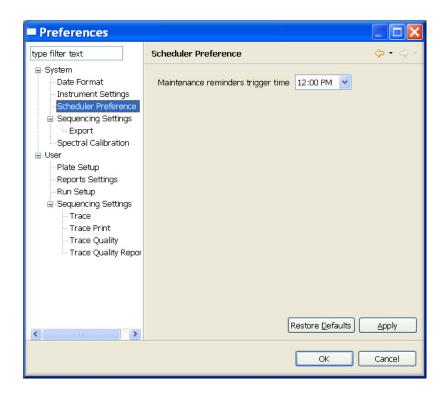
2. From the Left-hand pane, under Planned Maintenance, click Notifications Log

Click **(2)** on the top left-hand corner of the Notification Log for more information.

The Notification Log provides the following information on each event:

Notification	Description
Name	The name of the event.
Priority	The event priority.
Notification Date	The date of notification.
Status	The current status of the event.
User	The name of the user.
Acknowledge Date/Time	The date and time when the event was acknowledged.
Description	The description for the event.

Notification time is determined in the Preferences. From the Dashboard, click Preferences, to open the Preferences dialog box, click Scheduler Preference, and follow the prompts.



# Instrument operational procedures

The day-to-day operation of the instrument involves performing the following tasks.

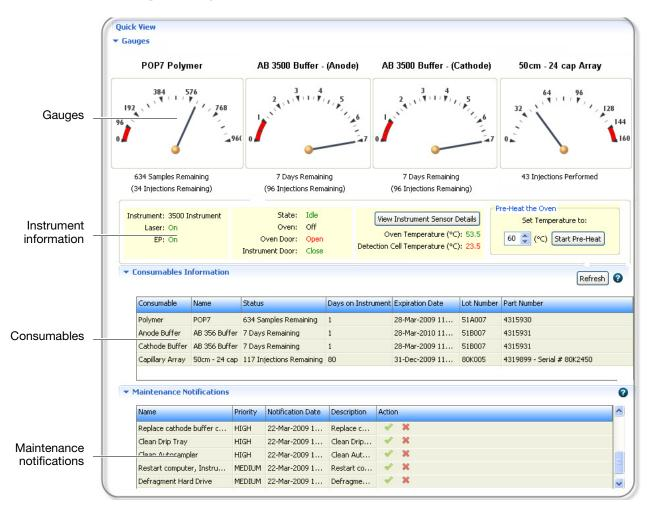
### Check consumables on the Dashboard

- Change the Anode Buffer Container (ABC)
- Change the Cathode Buffer Container (CBC)
- Change the polymer
- Use the Conditioning Reagent
- Fill Capillary Array with fresh polymer
- Remove bubbles

The Quick View section of the Dashboard provides the necessary information that you need to operate the instrument.

The information shown within the Quick View is generated automatically, via the Radio Frequency Identification (RFID) reader.

Use the information presented to you in the Quick View section before and after performing a maintenance task.



# Change the anode buffer container (ABC)

For the following hazard(s), see the complete safety alert descriptions in "Specific chemical alerts" on page 333.

# WARNING! CHEMICAL HAZARD. Anode Buffer Container (ABC).

For details see "Instrument reagents and consumables" on page 9.

Contamination might cause poor-quality data. To prevent the contamination, use genuine packaged polymer, anode buffer, cathode buffer and conditioning reagent.

- 1. Remove the ABC from storage.
- 2. Check for expiration date on the ABC label to make sure it is not expired prior to or during intended use.
- 3. Allow refrigerated ABC to equilibrate to ambient temperature prior to first use. Do not remove the seal until you have completed step 5, below.

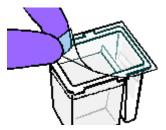
**IMPORTANT!** Ensure that all the buffer is moved to the larger side of the ABC prior to removing the seal.

4. Verify that buffer level is at or above the fill line and check that seal is intact.

**IMPORTANT!** Do not use if buffer level is too low or seal has been compromised. A fill tolerance of  $\pm 1$  mm is acceptable.

- Tilt the ABC slightly (as shown in the figure below) to make sure most of 1× buffer is in the larger side of the container. There should be less than 1 ml of 1× buffer remaining in the smaller side of the container.
- 6. Verify that the buffer is at the fill line.
- 7. Peel off the seal at the top of the ABC.





8. Place the ABC into the Anode end of the instrument, below the pump.

**IMPORTANT!** The RFID label must be facing the instrument (and not you) to ensure that the RFID information is read accurately by the instrument.



9. Close the instrument door to re-initialize.

**Note:** If you do not close the instrument door to re-initialize, you need to click **Refresh** from the Dashboard.

- 10. Click **Refresh** from the Dashboard to update the screen.
- 11. Check the Quick View section of the Dashboard for updated status after changing the ABC.

### Change the cathode buffer container (CBC)

For the following hazard(s), see the complete safety alert descriptions in "Specific chemical alerts" on page 333.

### WARNING! CHEMICAL HAZARD. Cathode Buffer Container (CBC).

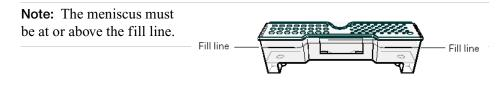
For details see "Instrument reagents and consumables" on page 9.

Contamination might cause poor-quality data. To prevent the contamination, use genuine packaged polymer, anode buffer, cathode buffer and conditioning reagent.

Use genuine parts and reagent. The use of inappropriate parts, or reagents, causes poor-quality data or damage the instrument.

- 1. Remove the CBC from storage.
- 2. Check for expiration date on the CBC label to make sure it is not expired prior to or during intended use.
- 3. Allow refrigerated CBC to equilibrate to ambient temperature before use.
- 4. Wipe away condensation on the CBC exterior with a lint-free lab cloth.
- 5. Verify that buffer level is at or above the fill line and check that seal is intact.

**IMPORTANT!** Do not use if buffer level is too low or seal has been compromised. A fill tolerance of  $\pm 0.5$ mm is acceptable.



6. Tilt the CBC back and forth gently and carefully to ensure that the buffer is evenly distributed across the top of the baffles.

**Note:** If you do not tilt the CBC back and forth, the buffer sticks to the baffles, due to surface tension.

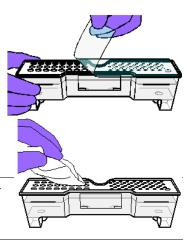
- 7. Verify that the buffer is at or above the fill line.
- 8. When ready to install CBC, place the container on a flat surface (such as a lab bench) and peel off the seal.
- 9. Wipe off any buffer on top of the CBC with a lint-free cloth. Ensure that the top of the container is dry.

**IMPORTANT!** Failure to perform this action may result in an arcing event and termination of the run.

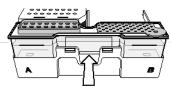
- 10. Place the appropriate septa on both sides of the CBC.
  - a. Align the buffer septa (the part that is symmetrical) over the 24 holes of the CBC.
  - b. Push the septa lightly into the holes to start and then push firmly to seat the septa.
- 11. Install the CBC on the autosampler.

**Note:** When properly installed, it will click on the autosampler as the tabs are snapped in place.

- 12. Close the instrument door to re-initialize.
- 13. Click **Refresh** from the Dashboard to update the screen.
- 14. Check the Quick View section of the Dashboard for updated status after changing the CBC.







### Check stored capillary arrays

**IMPORTANT!** Wear appropriate protection, including gloves, laboratory goggles, and coat whenever you work with the fluids used on this instrument, or parts that may come into contact with these fluids.

When the capillary array is installed, electrodes at the bottom are inserted on the CBC. The electrodes at the top connect with the polymer delivery pump. Applied Biosystems recommends you keep the electrodes on the bottom in the tray with  $1 \times$  running buffer. For details see "Instrument reagents and consumables" on page 9.

**IMPORTANT!** Keep the loading-end of the capillary array in  $1 \times$  running buffer to prevent the polymer from drying in the capillaries. If fluid level is low, add distilled water (DI) to buffer solution.

Refer to the Install capillary wizard for instructions on how to store the capillary array.



1X running buffer and distilled water (DI)

### Flush the water trap (pump trap)

The water trap must be flushed once per month to prolong the life of the pump and to clean any diluted polymer.

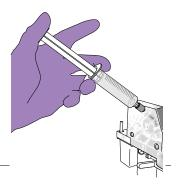
Flush with either distilled or deionized water and ensure that the water flows into the overflow container. Dispose the excess water (inside the overflow container). See "General chemical safety" on page 328.

Note: Leave the trap filled with either distilled or deionized water.

1. Fill the supplied 20 mL, all-plastic Luer lock syringe (in the PDP Cleaning kit, 4359572) with distilled or deionized water. Expel any bubbles from the syringe.

**IMPORTANT!** Do not use a syringe smaller than 20 mL. Doing so may generate excessive pressure within the trap.

- 2. Attach the syringe to the forward-facing Luer fitting at the top of the pump block. Hold the fitting with one hand while threading the syringe onto the fitting with the other hand.
- 3. Open the Luer fitting by grasping the body of the fitting and turning it to loosen. Attached syringe and turn counterclockwise approximately one-half turn.



IMPORTANT! DO NOT USE EXCESSIVE

FORCE when you push the syringe plunger as this may damage the trap seals. Take approximately 30 seconds to flush 5 mL of either distilled or deionized water through the trap.

**Note:** Because the water trap volume is approximately  $325 \ \mu$ L, a relatively small volume of water is adequate for complete flushing. However, a larger volume only improves flushing as long as force and flow rate are kept within the limits given above.

- 4. Remove the syringe from the Luer fitting. Hold the fitting with one hand while turning the syringe counterclockwise with the other hand.
- 5. Close the Luer fitting by lightly turning clockwise until the fitting seals against the block.

### Routine instrument cleaning

**IMPORTANT!** Wear appropriate protection, including gloves, laboratory goggles, and coat whenever you work with the fluids used on this instrument, or parts that may come into contact with these fluids.

- 1. Ensure the oven and instrument doors are closed.
- 2. Press the Tray button on the front of the instrument to move the autosampler to the forward position.

**IMPORTANT!** Use the cleaning agents as described in this manual, only. Use of cleaning agents not described in this manual can impair the instrument. Please contact your local Life Technologies sales office if you have any questions.

- 3. Wipe off any liquid on or around the autosampler using a lint-free tissue.
- 4. Clean off any polymer build-up crystals on the instrument, including the capillary tips, with deionized water and lint-free tissue.
- 5. Clean the array plug.
- 6. Clean out the drip trays with deionized water, or ethanol (absolute), and lint-free tissue.

Note: The drip tray can be removed.

### Move and level the instrument

**CAUTION! PHYSICAL INJURY HAZARD**. Do not attempt to lift the instrument or any other heavy objects unless you have received related training. Incorrect lifting can cause painful and sometimes permanent back injury. Use proper lifting techniques when lifting or moving the instrument. Two or three people are required to lift the instrument, depending upon instrument weight.

- 1. Remove the following components from the instrument:
  - Any plate assemblies from the autosampler.
  - CBC from the autosampler.
  - Capillary array: click **Shutdown the Instrument in the Maintenance** Wizards. See "To shutdown the instrument" on page 253.
  - Anode buffer reservoir.
- 2. Switch off the circuit breaker on the back of the instrument.
- 3. Disconnect the power cord and the Ethernet cable.

**IMPORTANT!** While moving the instrument, avoid any shock or vibration.

- 4. Move the instrument.
- 5. Turn the instrument legs to level the instrument.

To move the instrument corner	Turn the leg	
up	right (clockwise)	
down	left (counterclockwise)	

# Use the Maintenance Wizards to perform operations

# **About Maintenance Wizards**

To activate the Maintenance Wizards from the Dashboard, click **Maintain Instrument** toggle key.

The Maintenance Wizards feature of the Data Collection software allows you to perform operations necessary for sustaining the instrument.

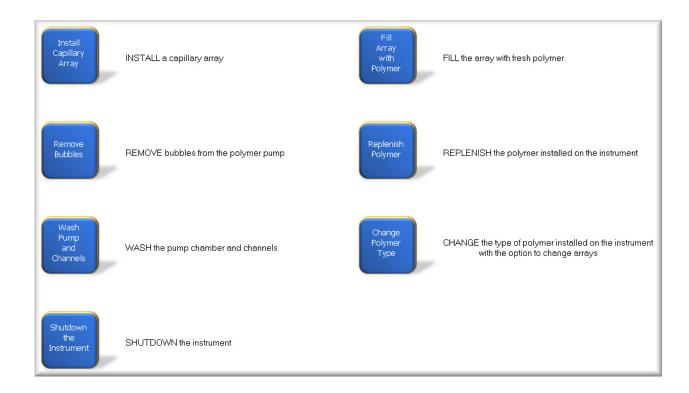
In no particular order, these operations include the following:

- Install a Capillary Array
- Remove bubbles from the polymer pump
- Wash the pump chamber and channels
- Fill the array with fresh polymer
- Replenish the polymer installed on the instrument
- Change the type of polymer installed on the instrument with the option to change the capillary array.
- Shutdown the Instrument.

**IMPORTANT!** Once started, Wizard operations cannot be canceled.

**IMPORTANT!** After performing a conditioning wash ensure that the buffer level inside the ABC is at or above fill line before proceeding to the next step except for the wash pump and channels wizard.





# **Replenish polymer**

**IMPORTANT!** Do not use a polymer pouch that has been installed on one type of instrument on another type of instrument. For example, if you install a new polymer pouch originally on a 3500 (8-capillary) instrument, do not subsequently use that same polymer pouch on a 3500xL (24-capillary) instrument, or vice versa. Doing so may result in a lower number of samples/injections than specified.

For the following hazard(s), see the complete safety alert descriptions in "Specific chemical alerts" on page 333.

**WARNING!** CHEMICAL HAZARD. POP- $4^{\text{TM}}$ , POP- $6^{\text{TM}}$ , and POP- $7^{\text{TM}}$  polymers.

For details see "Instrument reagents and consumables" on page 9.

If you are replacing the same polymer type only, follow the procedures below:

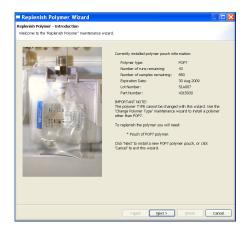
**IMPORTANT!** If you remove a polymer pouch for storage, place a Pouch Cap (PN 4412619) onto the pouch, then place an empty pouch (or conditioning reagent) on the connector to prevent desiccation of any residual polymer on the connector.

1. In the Maintenance Wizards screen, click Replenish Polymer.



Note: The Replenish Polymer Wizard takes 10 to 20 minutes to complete.

- 2. Follow the prompts in the Replenish Polymer Wizard window.
- 3. Click **Refresh** from the Dashboard to update the screen.
- 4. Check the Quick View section of the Dashboard for updated status after replenishing the polymer.



### Change polymer type

**IMPORTANT!** Do not use a polymer pouch that has been installed on one type of instrument on another type of instrument. For example, if you install a new polymer pouch originally on a 3500 (8-capillary) instrument, do not subsequently use that same polymer pouch on a 3500xL (24-capillary) instrument, or vice versa. Doing so may result in a lower number of samples/injections than specified.

For the following hazard(s), see the complete safety alert descriptions in "Specific chemical alerts" on page 333.



**IMPORTANT!** If you remove a polymer pouch for storage, place a Pouch Cap (PN 4412619) onto the pouch, then place an empty pouch (or conditioning reagent) on the connector to prevent desiccation of any residual polymer on the connector.

For details see "Instrument reagents and consumables" on page 9.

**IMPORTANT!** If the polymer dries on the fitment or in the pouch opening, the dried polymer prevents the pouch fitment from closing the internal cap properly. If that happens, the polymer pouch is no longer usable. When the pouch is removed, cover the fitment with a new, empty, or a conditioning pouch. To prevent drying, the pouch fitment must be covered with Pouch Cap (PN 4427991).

Note: Expired pouches cannot be used on the instrument.

- 1. Remove the polymer from storage 4 °C.
- 2. Allow refrigerated polymer to equilibrate to ambient temperature before use.
- 3. Check for expiration date on the pouch label to make sure it is not expired prior to use.

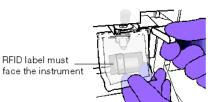
**IMPORTANT!** Do not use if the pouch and/or the label is damaged or the top seal is missing.

4. Peel off seal at the top of the pouch fitment.

**Note:** You may occasionally notice a tiny droplet of polymer inside the fitment (residual from the pouch filling process). This is **not** expected to cause any performance issues.

5. Slide the pouch fitment on to the slot of the lever assembly. Push the lever up to snap the pouch into the connector end of the instrument pump.

**Note:** The RFID label must be facing the instrument (and not you) to ensure that the RFID information is read accurately by the instrument.



- 6. If a partially used pouch is removed for later use, use the suggested cap to plug the fitment opening and store the pouch under recommended storage conditions.
- 7. From the Maintenance Wizards screen, click **Change Polymer Type**.



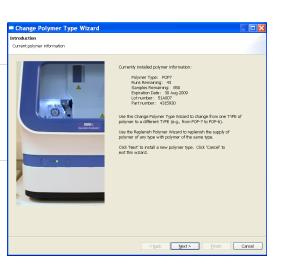
**IMPORTANT!** This feature allows you to change the type of polymer installed on the instrument with the option to change the Capillary Arrays.

Note: The Change Polymer Type Wizard takes 60 to 70 minutes to complete.

8. Follow the prompts in the Change Polymer Type Wizard window.

Note: Changing polymer requires the use of a Conditioning Reagent. See "Use the conditioning reagent" on page 250.

- 9. Click **Refresh** from the Dashboard to update the screen.
- 10. Check the Quick View section of the Dashboard for updated status after changing the polymer.



# Partially used polymer

**IMPORTANT!** Do not use a polymer pouch that has been installed on one type of instrument on another type of instrument. For example, if you install a new polymer pouch originally on a 3500 (8-capillary) instrument, do not subsequently use that same polymer pouch on a 3500xL (24-capillary) instrument, or vice versa. Doing so may result in a lower number of samples/injections than specified.

If a partially used pouch is removed for later use, use the suggested Pouch Cap to plug the fitment opening and store the pouch under recommended storage conditions. The Pouch Cap is sold separately (4412619).

If you remove a polymer pouch for storage, place a Pouch Cap (PN 4412619) onto the pouch, then place an empty pouch (or conditioning reagent) on the connector to prevent desiccation of any residual polymer on the connector. If the polymer dries on the fitment or in the pouch opening, the dried polymer prevents the pouch fitment from closing the internal cap properly. If that happens, the polymer pouch is no longer usable.

**IMPORTANT!** Follow the instructions in the wizard to ensure the proper installation and operation of the pouch and the instrument.

# Wash the pump chamber and channels

Note: The Wash Pump and Channels wizard takes over 40 minutes to complete.

1. From the Maintenance Wizards screen, click **Wash Pump** and Channels.



2. Follow the prompts in the Wash Wizard window.



### Use the conditioning reagent

For details see "Instrument reagents and consumables" on page 9.

**IMPORTANT!** Expired pouches cannot be used on the instrument. Once installed on the instrument, the pouch is good for a one-time use, only.

The use of the conditioning reagent is dictated by the instrument wizards.

Contamination might cause poor-quality data. To prevent the contamination, use genuine packaged polymer, anode buffer, cathode buffer and conditioning reagent.

Use genuine parts and reagent. The use of inappropriate parts, or reagents, causes poor-quality data or damage the instrument.

Refer to Chapter 3, Set Up and Run for instructions on priming the pump and initiating the run.

The Quick View section of the Dashboard provides the necessary information that you need for using the Conditioning Reagent.

Note: Install the pouch only when requested to do so by the wizard.

#### To place the conditioning reagent on the instrument

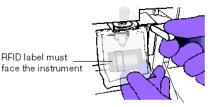
1. Check for expiration date on the label to make sure it is not expired prior to use.

**IMPORTANT!** Do not use if pouch/label is damaged or top seal is missing.

- 2. Peel off the seal at the top of the conditioning reagent pouch fitment.
- 3. Insert the pouch fitment on to the slot of the pump lever mechanism. Push the lever up to snap the pouch into the connector end of the instrument pump.

**Note:** The RFID label must be facing the instrument (and not you) to ensure that the RFID information is read accurately by the instrument.

- 4. Follow the wizard for further instructions.
- 5. Click **Refresh** from the Dashboard to update the screen.



 Check the Quick View section of the Dashboard for updated status after changing the Conditioning Reagent.

### Fill capillary array with fresh polymer

For the following hazard(s), see the complete safety alert descriptions in "Specific chemical alerts" on page 333.

# **WARNING!** CHEMICAL HAZARD. POP- $4^{\text{TM}}$ , POP- $6^{\text{TM}}$ , and POP- $7^{\text{TM}}$ polymers.

For details see "Instrument reagents and consumables" on page 9.

The filling of the capillary array with fresh polymer is dictated by the instrument wizards.

- 1. To fill capillary array with fresh polymer (same type of polymer), click **Fill the Array with fresh Polymer**.
  - er.



- 2. Follow the prompts in the Fill Array Wizard window.
- 3. Click **Refresh** from the Dashboard to update the screen.
- 4. Check the Quick View section of the Dashboard for updated status after filling of the Capillary Array with fresh polymer.



### Remove bubbles from the polymer pump

Remove bubbles from the polymer pump fluid path before each run. See "Daily instrument maintenance tasks" on page 230 for more information.

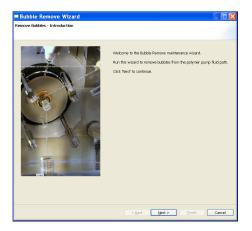
**IMPORTANT!** Wear gloves while handling polymer, the capillary array, septa, or CBC.

1. To remove bubbles from the polymer pump fluid path that travel from the polymer pouch through the pump, array port, and the Anode Buffer Container, click **Remove Bubbles**.

Remove Bubbles from the polymer pump

Note: The Bubble Remove Wizard takes 5 to 15 minutes to complete.

- 2. Follow the prompts in the Bubble Remove Wizard window.
- 3. Check the Quick View section of the Dashboard for updated status of the polymer pouch after removing bubbles from the polymer pump fluid path.



### To change the capillary array

**CAUTION! SHARP** The load-end of the capillary array has small but blunt ends and it could lead to piercing injury.

**IMPORTANT!** Check the loading-end header to ensure that the capillary tips are not crushed or damaged.

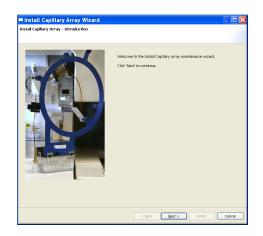
For details see "Instrument reagents and consumables" on page 9.

1. From the Maintenance Wizards screen, click **Install Capillary** Array.



Note: The Install Capillary Array Wizard takes 15 to 45 minutes to complete.

- 2. Follow the prompts in the Install Capillary Array Wizard window.
- 3. Check the Quick View section of the Dashboard for updated status of the capillary array.



# To shutdown the instrument

Use the Instrument Shutdown Wizard for short- and long-term shutdown.

1. From the Maintenance Wizards screen, click **Shutdown the Instrument**.

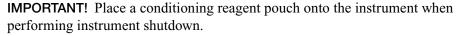
Note: The Instrument Shutdown Wizard takes 60 minutes to complete.

2. Follow the prompts in the Instrument Shutdown Wizard window.

Perform the appropriate shutdown procedure based on the information in the following table:

Instrument Shutdown Wizard	
Instrument Shutdown Wizard - Introduction	
() ····	The valued will help sporessore the internet for enderself protot of draws (grader the 3-coles), The optimizer area in more all in the proton in the optimizer area in the optimizer and the optimizer and the optimizer and area optimizer and the optimizer and area optimizer and area optimizer and area optimizer and area optimizer and area optimizer and area optimizer and area optimizer and area optimizer and
	<back mext=""> Einish Cancel</back>

SHUTTOWN



If the instrument will be unattended for	Perform this shutdown procedure
no more than 1 week	No action is required.
1 to 2 weeks	<b>IMPORTANT!</b> Keep the load-end of the capillary array in 1X buffer to prevent the polymer from drying in the capillaries. If fluid level is low, add DI water to buffer solution. Install the new CBC when ready to resume runs.
for more than 2 weeks	Long-term.
	See below for long-term instrument shutdown.

# Computer maintenance

This section lists the common tasks required to maintain the computer for your 3500 or 3500xL analyzer in good working condition.

For the computer troubleshooting issues, see Appendix E, "Troubleshoot" on page 299.

#### Uninstall the software

When you uninstall the software, you are prompted to back up the datastore (the directory that contains all library items you created, such as plates and protocols).

**IMPORTANT!** Do not back up the datastore to the installation directory. The installation directory is deleted during the uninstall.

#### Archive, purge, and restore data

- Archive Makes a copy of the data in an external file that you can save in another location.
- **Purge** Allows you to delete (purge) user-created items stored in the library. Factory-provided items are not purged. You have an option to archive the items, also.
- **Restore** Restores archived data back to the system.

**IMPORTANT!** These functions affect items stored in the library (datastore). These functions do not affect sample data files.

- **Frequency** Applied Biosystems recommends that you purge the library objects once every three months.
- Archive library This function archives items stored in the library. To archive audit records, see "Archive, purge, and restore audit records" on page 214.
  - 1. Access the Archive screen.



2. Specify the date category and range, then click OK.

	Date <ul> <li>Created</li> <li>Modified</li> <li>All</li> </ul>	
	Date Range from //  to //	
	OK Cancel	
	3. Specify a location and file name for the archive (.dsz) file, then click Save.	
	A message is displayed when the archive is complete.	
	<b>IMPORTANT!</b> Do not specify x:\Applied Biosystems\3500\datastore as the archive location. If you do so, your archive can be deleted if you uninstall the software and do not back up the datastore.	
	If you specify a location to which you do not have permission to save, a warning message is displayed and gives you the option to save in another location.	ng
Archive data files	There are two ways to archive the data files	
	1. Start > Control Panel > System and Maintenance > Backup and Restore Center	er

Start 
 Control Panel 
 System and Maintenance 
 Backup and Restore Center
 OR

Programs > Accessories > System tools > Backup

2. Use either Back up File folder or Back up Computer options.

**Note:** If you export audit records for samples that are not in their original location (samples have been deleted or moved), an error message is displayed. Return sample data files to their original location, then export again.

- **Restore** This function restores items stored in the library. To restore audit records, see "Archive, purge, and restore audit records" on page 214.
  - 1. Access the Restore function.



2. Select the archive (.dsz) file to restore, then click **Open**.

If the archive file contains items that exist in the system, a message is displayed.



3. Select an option to continue.

A message is displayed when the restore is complete.

- **Purge** This function purges (deletes) items stored in the library. To purge audit records, see "Archive, purge, and restore audit records" on page 214.
  - 1. Access the Purge function.



- 2. Click **Yes** in the Purge warning message stating that you are about to permanently delete all files in the library.
- 3. Specify the date category and range, then click OK.
- 4. Click Yes in the Purge warning message.

A message is displayed when all records are deleted.

### Monitor disk space

Ensure that you have sufficient drive space by regularly:

- Archiving data
- Deleting unneeded files
- Emptying the trash
- Defragmenting the drives

Hard disk and<br/>statusManually check available disk space on Drive D.To check the status, go to My Computer ▶ right-mouse click on C drive ▶ Select<br/>Properties ▶ Click General tab.

		Local Disk (C:) Properties		
	Note: The Data Collection software will prompt you when it is	General Tool: Hardware Sharing Security Quota		
	70-75% full. At 78% full, the software will not start a run.	Type: Local Disk File system: NTFS		
		Used space: 31,548,768,256 bytes 23,3 GB Free space: 48,393,825,408 bytes 45,0 GB Cepority: 73,908,593,694 bytes 74,4 GB		
	If there is insufficient space:			
	• Archive the sample files.	Drive C Disk Dearup		
	• Delete the sample file data from the drive D and empty the contents of the Recycle Bin.	Compress diver to save dui space Calculate indexing Service to yolds the duit for fast the searching OK Cancel Archive		
Defragment the computer hard drive	This option can be set as a reminder in the scheduler. The fragmer decreases the performance of both the Data Collection software an operating system. Programs take a longer time to access files by per- search operations of the fragments.	nd the computer		
	Go to Start > Programs > Accessories > System Tools > Disk Defragmenter and follow the prompts.			
	Note: You can click Analyze to see if you should defragment or r	lot.		
Check available space on all drives	Before a run, the Data Collection software checks free disk space. disk space is not available to store the data, the Data Collection so following message:	•		
	Remove data: the drive is getting full			
	View the errors that appear for generated errors and in the Event I Appendix E, "Troubleshoot" on page 299.	log window. See		
	Also, check the status light in the bottom left-hand corner of the d window to see if it flashes red.	ata collection		

# **Review the Maintenance Notifications Log**

The Notifications Log is a history of all notifications messages and the action taken for the task (completed or dismissed). You can use this option to review a previous run information.

The Dashboard provides you with a list of current routine and maintenance notifications, as explained below.

Multi-column sorting is supported (see "Multi-column sorting" on page 72).

To go to the Notifications Log from the Dashboard:

#### 1. Click Maintain Instrument

2. From the Left-hand pane, under Planned Maintenance, click Notifications Log

Click ② on the top left-hand corner of the Notification Log for more information.

Notification	Description
Name	The name of the event.
Priority	The event priority.
Notification Date	The date of notification.
Status	The current status of the event.
User	The name of the user.
Acknowledge Date/Time	The date and time when the event was acknowledged.
Description	The description for the event.

The Notification Log provides the following information on each event:

Notification time is determined in the Preferences. From the Dashboard, click Preferences, to open the Preferences dialog box, click Scheduler Preference, and follow the prompts.

# Service Log

The Service Log is a record of instrument service, and it is used and completed by the Applied Biosystems service engineer at the time of service.

To go to the Service Log from the Dashboard:

#### 1. Click Maintain Instrument

2. From the Left-hand pane, under Planned Maintenance, click Service Log

Click **(2)** on the top left-hand corner of the Service Log for more information.

The Service Log screen contains a history of all the service events that have occurred on the system, starting with the most recent event, and provides the following information on each event:

Event	Description
Ticket Number	The number assigned to the event.
Service Type	The type of service requested.
Event Occur Date	The date that the event took place.
Service Start Date	The date that the service started.
Service End Date	The date that the service ended.
Service Engineer	The name of the service engineer.
Reason	The reason for logging the event.
Comments	Any additional comments.

# Sequencing analysis reagents

**Note:** For more details see the product insert included in the product package.

The following table shows all the reagents for sequencing analysis.

#### Table 27 Sequencing analysis reagents

Name	Part Number	Storage Conditions	On-instrument Shelf-life at Environmental Temperature
BigDye <sup>®</sup> Terminator (BDT) v3.1 Cycle Sequencing Kit 24 reactions	4337454	-15 °C to -25 °C	24 hours
BigDye <sup>®</sup> Terminator (BDT) v3.1 Cycle Sequencing Kit 100 reactions	4337455	-15 °C to -25 °C	24 hours
BigDye <sup>®</sup> Terminator (BDT) v3.1 Cycle Sequencing Kit 1000 reactions	4337456	-15 °C to -25 °C	24 hours
BigDye <sup>®</sup> Terminator (BDT) v3.1 Cycle Sequencing Kit 5000 reactions	4337457	-15 °C to -25 °C	24 hours
BigDye <sup>®</sup> Terminator (BDT) v1.1 Cycle Sequencing Kit 24 reactions	4337449	-15 °C to -25 °C	24 hours
BigDye <sup>®</sup> Terminator (BDT) v1.1 Cycle Sequencing Kit 100 reactions	4337450	-15 °C to -25 °C	24 hours
BigDye <sup>®</sup> Terminator (BDT) v1.1 Cycle Sequencing Kit 1000 reactions	4337451	-15 °C to -25 °C	24 hours
BigDye <sup>®</sup> Terminator (BDT) v1.1 Cycle Sequencing Kit 5000 reactions	4337452	-15 °C to -25 °C	24 hours

#### Table 28Sequencing standards

Name	Part Number	Storage Conditions	On-instrument Shelf-life at Environmental Temperature
BigDye® Terminator (BDT) v3.1 Sequencing Standard (long read)	4404312	-15 °C to -25 °C	24 hours

#### Table 28 Sequencing standards

Name	Part Number	Storage Conditions	On-instrument Shelf-life at Environmental Temperature
BigDye® Terminator (BDT) v1.1 Sequencing Standard (long read)	4404314	-15 °C to -25 °C	24 hours
BigDye <sup>®</sup> Terminator (BDT) v3.1 Matrix Standard	4336974	2 °C to 8 °C	24 hours
BigDye <sup>®</sup> Terminator (BDT) v1.1 Matrix Standard	4336824	2 °C to 8 °C	24 hours

# Fragment and HID analysis reagents

**Note:** For reagent or consumable shelf-life expiration date, see the package label.

The following table shows all the reagents for fragment and HID analysis.

Table 29 Fragment analysis HID standards

Name	Part Number	Storage Conditions	On-instrument Shelf-life at Environmental Temperature
Fragment Analysis Matrix Standards (5-Dye) -DS-02	4323014	2 to 8°C	24 hours
Fragment Analysis Matrix Standards (4-dye) - DS-32	4345831	2 to 8°C	24 hours
Fragment Analysis Matrix Standards (5-Dye) -DS-33	4345833	2 to 8°C	24 hours
Fragment Analysis Installation kit (5-Dye) -DS-33	4376911	2 to 8°C	24 hours
GS120LIZ Size Standard	4322362	2 to 8°C	24 hours
GS500ROX Size Standard	401734	2 to 8°C	24 hours
GS600 LIZ Size Standard v2 (for Normalization)	4408399	2 to 8°C	24 hours
GS1200 LIZ Size Standard	4379950	2 to 8°C	24 hours

# Sequencing analysis dye sets for all applications

Note: For reagent or consumable shelf-life expiration date, see the package label.

The following table shows all the dye sets for various applications.

Table 30 Dye Sets for various applications

Dye Set	Application Name
E (v1.1 BigDye <sup>®</sup> Terminator)	Rapid DNA sequencing
Z (3.1 BigDye® Terminator)	DNA sequencing

# Fragment analysis dye sets for all applications

Note: For reagent or consumable shelf-life expiration date, see the package label.

The following table shows all the dye sets for fragment analysis.

Table 31Fragment analysis dye sets

Dye Set	Application
E5	SNaPshot <sup>®</sup> kit
G5	DNA sizing for 5-dye chemistry
J6	DNA sizing for 6-dye chemistry
F	DNA sizing for 4-dye chemistry
Any dye	DNA sizing

# HID analysis dye sets

AmpF <b>/STR<sup>®</sup> Kits</b>	Dye set (use with HID Fragment Analysis 36_POP4 run module)
<ul> <li>4-dye:</li> <li>COfiler<sup>®</sup></li> <li>Profiler Plus<sup>®</sup></li> <li>Profiler Plus<sup>®</sup> <i>ID</i></li> <li>SGM Plus<sup>®</sup></li> <li>Other 4-dye kits</li> </ul>	F
5-dye: Identifiler <sup>®</sup> Minifiler <sup>™</sup> SEfiler <sup>™</sup> Plus SinoFiler <sup>™</sup> Yfiler <sup>®</sup> Other 5-dye kits	G5

#### Table 32 AmpF/STR Kit Table

# **Run modules**

# Capillary array and polymer (sequencing analysis run modules)

Decide what combination of capillary array and polymer matches your resolution and performance specifications, from the table below.

Table 33 Capillary array and polymer (sequencing analysis run modules)

Run Module Type &	Config	uration	23 hours Throughput <sup>‡</sup>			Performance
Run Module Name	Capillary Length (cm)	Polymer Type	Run Time (min)	3500	3500xL	Contiguous Read Length (CRL) <sup>§</sup>
Rapid sequencing RapidSeq50_POP7	50	POP-7 <sup>™</sup>	≤40	≥280	≥840	≥500
Standard sequencing StdSeq50_POP6	50	POP-6 <sup>™</sup>	≤135	≥80	≥240	≥600
Fast sequencing FastSeq50_POP7	50	POP-7 <sup>™</sup>	≤65	≥168	≥504	≥700
Standard sequencing StdSeq50_POP7	50	POP-7 <sup>™</sup>	≤125	≥88	≥264	≥850
Short read sequencing ShortReadSeqPOP7	50	POP-7 <sup>™</sup>	≤30	≥368	≥1104	≥300
Rapid sequencing BigDye <sup>®</sup> XTerminator <sup>™</sup> RapidSeq_BDX_50_POP7	50	POP-7 <sup>™</sup>	≤40	≥280	≥840	≥500
Standard sequencing BigDye <sup>®</sup> XTerminator <sup>™</sup> StdSeq_BDX_50_POP6	50	POP-6 <sup>™</sup>	≤140	≥80	≥240	≥600
Fast sequencing BigDye <sup>®</sup> XTerminator <sup>™</sup> FastSeq_BDX_50_POP7	50	POP-7 <sup>™</sup>	≤65	≥168	≥504	≥700
Standard sequencing BigDye <sup>®</sup> XTerminator <sup>™</sup> StdSeq_BDX_50_POP7	50	POP-7 <sup>™</sup>	≤125	≥88	≥264	≥850
Short read sequencing BigDye <sup>®</sup> XTerminator <sup>™</sup> ShortReadSeq_BDX_POP7	50	POP-7 <sup>™</sup>	≤30	≥368	≥1104	≥300
Microbial Sequencing MicroSeq_POP7	50	POP-7™	≤125	≥88	≥264	≥850

Run Module Type &	Configu	uration	23 ho	Performance		
Run Module Name	Capillary Length (cm)	Polymer Type	Run Time (min)	3500	3500xL	Contiguous Read Length (CRL) <sup>§</sup>
Microbial Sequencing MicroSeq_POP6	50	POP-6 <sup>™</sup>	≤135	≥80	≥240	≥600

#### Table 33 Capillary array and polymer (sequencing analysis run modules) (continued)

‡ Throughput (Samples / Day): The total number of samples run in 23 hours (0.5 hour for User interaction and 0.5 hour for warm-up time).

§ The maximum number of contiguous bases in the analyzed sequence with an average QV ≥20, calculated over a sliding window 20 base pairs wide from an AB Long Read Standard sequencing sample. This calculation starts with base number 1. The read length is counted from the middle base of the 1st good window to the middle base of the last good window, where a "good" window is one in which the average QV ≥20.

# Capillary array and polymer (fragment and HID analysis run modules)

	Configuration		23 hours Throughput <sup>‡</sup>			Performance			
Run Modules Type & Run Modules Name	Capillary		Run Time (min)	3500	3500xL	Range <sup>§</sup>	Sizing Precision <sup>#</sup>		
	Length (cm)	Polymer Type					50bp- 400bp	401bp- 600bp	601bp- 1200bp
Fragment analysis FragmentAnalysis50_POP7	50	POP-7 <sup>™</sup>	≤40	≥280	≥840	≤40 to ≥520	<0.15	<0.30	NA <sup>‡‡</sup>
Fragment analysis FragmentAnalysis50_POP6	50	POP-6 <sup>™</sup>	≤100	≥112	≥336	≤20 to ≥550	<0.15	<0.30	NA <sup>‡‡</sup>
Long fragment analysis LongFragAnalysis50_POP7	50	POP-7 <sup>™</sup>	≤125	≥88	≥360	≤40 to ≥700	<0.15	<0.30	<0.45
HID HID36_POP4	36	POP-4 <sup>™</sup>	≤35	≥312	≥936	≤60 to ≥400	<0.15	NA <sup>‡‡</sup>	NA <sup>‡‡</sup>
HID HID36_POP7	36	POP-7 <sup>™</sup>	≤26	≥424	≥1272	≤60 to ≥400	<0.15	NA <sup>‡‡</sup>	NA <sup>‡‡</sup>
SNaPshot <sup>®</sup> SNaPshot50_POP7	50	POP-7™	≤30	≥376	≥1104	≤40 to ≥120	<0.50	NA <sup>‡‡</sup>	NA <sup>‡‡</sup>

#### Table 34 Capillary array and polymer (fragment and HID analysis run modules)

‡ Throughput (Samples / Day): The total number of samples run in 23 hours (0.5 hour for User interaction and 0.5hr for warm-up time).

- § Resolution Range: The range of bases over which the resolution (peak spacing interval divided by the peak width at halfmax in a GS600 or GS1200 LIZ size standard sample sized with a third order fit) is ≥1. The table shows the resolution range in ≥90% of samples.
- # Sizing Precision: Standard deviation of sizes for one allele in the DS-33 install standard sized with the GS600 LIZ size standard across multiple capillaries in the same run. For one injection to pass, 100% of the alleles in that injection must meet the intra-run sizing precision specifications. The table shows the sizing precision of 100% of alleles in ≥90% of samples.
- **‡**‡Not applicable because of the size of the fragments collected in the run.

# Perform secondary analysis on sequencing experiments

The Applied Biosystems 3500/3500xL Genetic Analyzers and 3500 Series Data Collection Software provide integration between the instrument and secondary sequencing analysis software applications— specifically SeqScape<sup>®</sup> Software v2.7 and MicroSeq<sup>®</sup> ID Software v2.2. Using auto-analysis, samples are loaded, sequencing data is generated, and basecalling along with secondary analysis is performed according to the protocols assigned to the plates prior to the run.

Software	Purpose
SeqScape®	A comprehensive resequencing tool designed to detect SNPs, profile mutations, perform medical sequencing, identify haplotypes, subtype pathogens, and confirm clone constructs.
MicroSeq <sup>®</sup> ID	A comparative sequencing tool for microbial identification of bacteria and fungi.

# Auto-analyze projects in the sequencing analysis software

Auto-analysis can only be performed on the same computer that collects the sample files, therefore SeqScape<sup>®</sup> or MicroSeq<sup>®</sup> ID Software must be co-installed and configured with the 3500 Series Data Collection Software on a Windows Vista<sup>®</sup> operating system. Automated basecalling occurs with KB<sup>TM</sup> Basecaller v1.4.1 (calls pure or mixed bases with quality values) and secondary analysis occurs with SeqScape<sup>®</sup> or MicroSeq<sup>®</sup> ID Software.

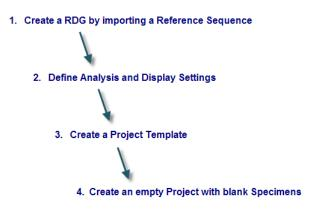
This procedure initially describes how to set up panels and bin sets in SeqScape<sup>®</sup> and then describes how to auto-analyze samples using the 3500 Series Data Collection Software. Once a run is complete, your data is seamlessly transferred into SeqScape<sup>®</sup> for analyzing, processing and reporting.

**Note:** For detailed information on setting up a MicroSeq<sup>®</sup> ID project to auto-analyze in the 3500 Series Data Collection Software, see the *MicroSeq<sup>®</sup> ID v2.2 Getting Started Guide*.

# Set up an auto-analysis project in SeqScape®

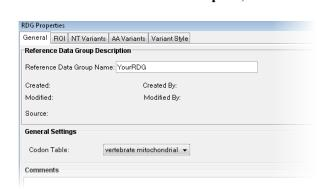
**IMPORTANT!** When using SeqScape<sup>®</sup> Software to auto-analyze results data from the 3500/3500xL analyzer, you must have v2.7 installed on the *same* computer as the 3500 Series Data Collection Software.

Set up a project in the secondary analysis software before starting a run on the 3500/3500xL analyzer. All analysis in SeqScape<sup>®</sup> occurs in a project. Create a project by following these steps:



Import a reference

Start the SeqScape<sup>®</sup> Software (<sup>™</sup><sub>seqScape</sub>), then select Tools → SeqScape Manager.
 Select the Reference Data Group tab, then click New and enter a name.



3. Select the **ROI** (Regions of Interest) tab, then click **Add Ref. Segment**.



- 4. Select the file you want to use as your reference file, then click Import.
- 5. In the **NT Variants** tab (RDG Properties), select the NT Variants that you want to add to the reference sequence, then click **Import** to import a tab-delimited variants file or a multi-aligned sequence (.fsta) file.

Note: When importing an amino acid variants file, use a tab-delimited format.

6. Click OK.

#### **Define settings** 1. Open **Tools** > **SeqScape Manager**.

- 2. Select the Analysis Protocols tab, then click New to enter a name.
- 3. Select the **Basecalling** tab, then select your Basecaller and Dye/Primer files.

Analysis P	rotocol Editor			
General	Basecalling	Mixed Bases	Clear Range	Filter
Baseca	-	271 B ben		_
DyeSet	/ Primer :			_
DT373	0POP7 (BDv3)	.mob		•
Proces	sed Data			
True	Profile			-
<ul> <li>Flat</li> </ul>	Profile		<b>Willian</b> a kinta	

**Note:** Unless a project requires a custom setting, keep the Processed Data, Ending Base and Quality Threshold settings at their default values.

- 4. In the Mixed Bases tab, specify the secondary analysis peak threshold for mixed base identification.
- 5. Keep the default settings for the other parameters listed in the Clear Range and Filter tabs, then click **OK**.
- 6. From the SeqScape Manager, select the **Analysis Defaults** tab, then click **New** and enter an Analysis Defaults Name.
- 7. Go to the **Sample** tab and select the Analysis Protocol you just created in the drop-down list.
- 8. Keep the default settings in the Project and Specimen tabs, then click Save. In most cases, you will want to keep the default Display Settings and continue with creating a project template in the SeqScape<sup>®</sup> Software.

Create a project	1. Open Tools > SeqScape Manager.
template	2. Select the <b>Project Templates</b> tab, then click <b>New</b> and enter a name for the template.
	3. In the Reference Data Group and Analysis Defaults drop-down lists, select the RDG and Analysis Default that you previously created.
	4. Keep the default Display Settings, then click <b>OK</b> .
	With the project template created, continue with adding your sample files.
Create an empty project	1. In the SeqScape <sup>®</sup> Software, select <b>File → New Project</b> .
	2. Name the project.
	3. In the Project Template list, select the project template that you previously created.
	4. When the project opens, click Add Specimen (Tools → New Specimen) to create as many blank specimens as you have in your project, then click OK.
	5. Close the SeqScape <sup>®</sup> Software.
	You are now ready to set up a run in the 3500 Series Data Collection Software specifying a SeqScape Protocol as your secondary analysis method.

# Set up a SeqScape plate in the 3500 Series Data Collection Software

Start the 3500 Series Data Collection Software

- 1. Start the Auto-Analysis Manager before starting the 3500 Series Data Collection Software.
- 2. Start the 3500 Series Data Collection Software ➤ Dashboard ➤



- 3. Name your new plate.
- 4. Select the Number of Wells, Plate Type as Sequencing, Capillary Length and Polymer associated with this plate for the current run.
- 5. (Optional) Enter your name as Owner, a barcode and description for the plate.

Specify autoanalysis for secondary analysis 1. Check **Perform Auto-Analysis** (right side of the Plate Details section), to expand the Secondary Analysis section.

📮 3500 Data Collection Software			
Dashboard Edit 🔻	Librar	ry Maintenance Tools 🔻 M	lanage 🔻 Preferences Help 🔻 Log Ou
🧱 Plate Name:	📖 New Plate 🔻 📄 Open Plate 🛛 🕞 Save Plate 🗸 📰 Close Plate	属 Start Run	
AB Applied Biosystems	Plate Details		0
Setup			
Define Plate Properties	* Name: New SeqScape Run	Owner:	
Assign Plate Contents	* Number of Wells:   96 96-FastTube 384	Barcode:	
Run Instrument Load Plates for Run Preview Run Monitor Run	* Plate Type: Sequencing • * Capillary Length: 50 • cm * Polymer: POP7 •	Description:	first sequence on 3500 🔹
Review Results View Sequencing Results View FragmenthID Results	Secondary Analysis      Software Type: SeqScape     Software Location: SeqScape_3500-FVTEST     Username:         *Password:      Assign Plate	e Contents	Perform Auto-Analysis 🛛

2. Confirm **SeqScape** auto-populates as the Software Type.

**Note:** If SeqScape does *not* appear in the drop-down list under Software Type, check your installation. Secondary analysis software must be installed correctly before the 3500 Series Data Collection Software is automatically listed as a selection.

3. Confirm Your computer name auto-populates as the Software Location.

**IMPORTANT!** For auto-analysis to be successful, the secondary analysis protocol must match the software location set here.

- 4. Enter your Username and Password for auto-analysis access to the secondary analysis software.
- 5. Click Save Plate > Save to save your plate with these settings, then Assign Plate Contents to advance to the next screen.

#### **Assign plate contents** When assigning plate contents, you are assigning assays, file name conventions and results group to be associated with your auto-analysis.

Set up an assay 1. In the Assign Plate Contents screen, go to the Assays box and select either Create New Assay or Add From Library.

Assays	
	Actions 🔻
Add from Library	
Create New Assay	
create rear rosay	

2. Name your assay in the Setup an Assay dialog box.

Note: (Optional) Select a color for this assay to display with in the Plate View.

3. Select an Instrument Protocol to apply to the assay.

**Note:** For more instruction on setting up an instrument protocol, see "Create a new instrument protocol" on page 165.

4. Select a Basecalling Protocol to apply to the assay.

**IMPORTANT!** Make sure your basecalling settings match the Analysis Settings specified in SeqScape.

**Note:** For more instruction on setting up a Basecalling protocol, see "Create a new basecalling protocol" on page 174.

5. Create a new sequencing analysis protocol to apply to the samples by clicking **Create New**.

📜 Create New Sequ	encing Analysis Proto	col		<b>X</b>
	ncing Analysis Pr			4
😢 Protocol Name	"Test" already exists in	the Library.		The second secon
				0
				-
* Protocol Name:	Test		Cocked	
Description:				
	Application Type:	Sequencing	-	
* Seconda	ary Analysis Software:	SeqScape	w	
* Secondary Analy	sis Software Instance:	SeqScape_fospradhaaaD03	•	
Properties				
•	* Project:	p53 v2	•	
	* Project Template:	pos_exon/	•	
	* Specimen:	12586	•	
Close				Save

6. Name your new sequencing analysis protocol, then select your specimens one by one, clicking **Save** after each specimen.

**IMPORTANT!** Each SeqScape protocol has one specimen, so you will need to create multiple protocols for multiple specimens. If you have multiple protocols, you will have multiple assays, as each assay is associated with one secondary analysis protocol.

**Note:** For more instruction on setting up a secondary analysis protocol, see "Create a new sequencing analysis protocol" on page 189.

New SeqScape Run - (	create riew Assay				-
etup an Assay					E
Assay Name: New	SeqScape Assay		Locked	Color:	Dark Cyan
pplication Type: Sequ	encing 👻				
Protocols					
	multiple instrume	nt protocols to this assay? 🔘 No	Yes		
		nt protocols to this assay?  ON DxFastSeq50_POP7_1	🔘 Yes	▼ Edit	Create New
Do you wish to assign	:			<ul><li>▼ Edit</li><li>▼ Edit</li></ul>	
Do you wish to assign * Instrument Protoco	l:	BDxFastSeq50_POP7_1			
Do you wish to assign * Instrument Protoco * Basecalling Protoco	l:	BDxFastSeq50_POP7_1 3500-MDC-Basecaling Prot		▼ Edit	Create New

7. Click Apply to Plate, then Save to Library if you want to use this assay again.

#### 8. Click Close.

AB Biosystems		View 🧱 Ta											
Setup										-			
Define Plate Properties			2	Show In W	ells 🔻 🔚	Select Wells	• An	ay Selection	Row	<table-of-contents> Column</table-of-contents>	200m	in 🔝 Zoom	Out 💽 Fit
Assign Plate Contents		1	2	3	4	5	6	7	8	9	10	11	12
Run Instrument	А	S sample100 A01	S sample108 A02	S sample116 A03	S sample124 A04	S sample132 A05	S sample140 A06	S sample148 A07	S sample156 A08	S sample164 A09	S sample172 A10	S sample180 A11	S sample188 A12
Load Plates for Run Preview Run	В	S sample101 B01	S sample109 B02	S sample117 B03	S sample125 B04	S sample133 B05	S sample141 B06	S sample149 B07	S sample157 B08	S sample165 B09	S sample173 B10	S sample181 B11	S sample189 B12
Monitor Run	с	S sample102 C01	S sample110 C02	S sample118 C03	S sample126 C04	S sample134 C05	S sample142 C06	S sample150 C07	S sample158 C08	S sample166 C09	S sample174 C10	S sample182 C11	S sample190 C12
Review Results	D	S sample103 D01	S sample111 D02	S sample119 D03	S sample127 D04	S sample135 D05	S sample143 D06	S sample151 D07	S sample159 D08	S sample167 D09	S sample175 D10	S sample183 D11	s sample191 D12
View Sequencing Results	E	S sample104 E01	S sample112 E02	S sample120 E03	S sample128 E04	S sample136 E05	S sample144 E06	S sample152 E07	S sample160 E08	S sample168 E09	S sample176 E10	S sample184 E11	S sample192 E12
View Fragment/HID Results	F	S sample105 F01	S sample113 F02	C	S	C	S sample145 F06	S sample153 F07	C	S	C	S sample185 F11	S sample193 F12
	G	S sample106 G01	S sample114 G02	S sample122 G03	S sample130 G04	S sample138 G05	S sample146 G06	S sample154 G07	S sample162 G08	S	S sample178 G10	S sample186 G11	S sample194 G12
and the	н	S sample107	S sample115	S sample123	S sample131	S sample139	S sample147	S sample155	S sample163	S sample171	S sample179	S sample187	S sample195

9. Name your samples by highlighting the number of wells in your plate and naming the sample in Customize Sample Info box.

			+ Custo	mize Sample Info	
		•	Property	Value	
Assays	File Name Conver	tions	1: Sample		
Actions 🔻		Actions	Sample Name	sample100	
🗹 🙉 BDx Fast Seq Assay 📝 🗙			Sample Type	Sample	=
🖉 🖉 BDX Fast Seq Assay 🖉 🔼	As General File Name	2	2: Custom		
			User Defined Fi	•	
			User Defined Fi	e	
			User Defined Fi	e la	
			User Defined Fi	¢	
			User Defined Fi	e	-
		E.	•		•

**Note:** For more information on naming samples, see "Name samples in the Plate View" on page 70.

Specify FNC and RG 1. Specify a File Name Convention (FNC) and a Results Group (RG) to associate with your project.

**Note:** You can create a FNC with the specimen name as a part of your sample file name.

- 2. Highlight the wells of your plate configuration (Plate View) and check the box next to the appropriate FNC to apply it to your project.
- 3. Repeat for the Results Group.

**Note:** For more information on setting up a FNC see "Create a new file name convention" on page 151. For more information on setting up a RG, see "Create a new results group" on page 156.

4. Click ■ Save Plate ► Save.

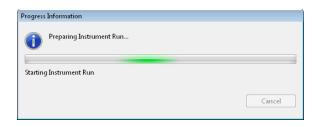
Plate View 🕎 Table View		
💽 Show In Wells 👻 🖃	Select Wells 🔻 📗 Array Selection 🛛 💽 Row 🚍 Co	lumn 🔢 Zoom In 🔝 Zoom Out 💽 Fit 💡
	5 6 7 8	
		9 10 11 12 5 5 5 5 5 5
	F \$ \$ \$ \$ \$ \$ \$ \$ \$ \$ \$ \$ \$ \$ \$ \$ \$ \$ \$	
	રક્રિક્સક્સક્સ્ક્સ્ક્સ્ટ્સ્ક્સ્ટ્સ્ક્સ્ટ્સ્ટ્	
	ይአቀኝኛ ላጅ ናይአቀኝያ ላጅ ናይአቀኝያ ላጅ ናይቀኝያ ላጅ ና	
	LS &A &&&axax& & & &&&&&&&&&&&&&&&&&&&&&&&	
	LS && 4550x45 && 4550x45 && 4550x45 && 4550x45	ት ትምምምምምምምምምምምምምምምምምምምምምም
F &A Assay & S &A Assay & S &A Assay & S &A Assay & S	LS &A 4550X &S &A 4550X &S &A 4550X &S &A 4550X &S	**************************************
G &A Assay & S &A Assay & S &A Assay & S &A Assay &	S & Assaves & Assaves & Assaves & Assaves & Assaves	A Assave & A Assave & A Assave & A Assave
H & Assay, A Si & Assay, A Si & Assay, A Si & Assay, A Si	Si 🗛 Assay, 4 Si 🗛 Assay, 4 Si 🗛 Assay, 4 Si 🗛 Assay, 4 Si	🗛 Assay,4 Si 🗛 Assay,4 Si 🗛 Assay,4 Si 🍇 Assay,4 🤟
S Sequencing		
s sequencing		
Name: AA Seq Plate 4 SeqScape	Barcode: sqa	
		0
Assays	File Name Conventions	Results Groups
Actions 🔻	Actions 🔻	Actions
● 🖌 AA Assay 4 SeqScape 📝 🗙	💿 🗹 TEST FNC 🛛 📝 🛛 🗙	● 🗹 TEST_RG 🛛 📝 🗵
AA Assay 4 SeqScape 2 X		● 🗹 TEST_RG 🛛 🖉 🔀
	● 🗹 TEST FNC 📝 🗙	● 🗹 TEST_RG 🥢 🔀
<ul> <li>AA Assay 4 SeqScape</li> <li>AA Assay 4 SeqScape</li> </ul>	● Z TEST FNC Z	Action:
		● ☑ TEST_RG 🔗 🔀
		● ☑ TEST_RG

- 5. Click Link Plate for Run.
- 6. Click **Create Injection List**, then click **OK** after the instrument performs its validations.

#### Start the autoanalysis run

Click Start Run to begin your auto-analysis.

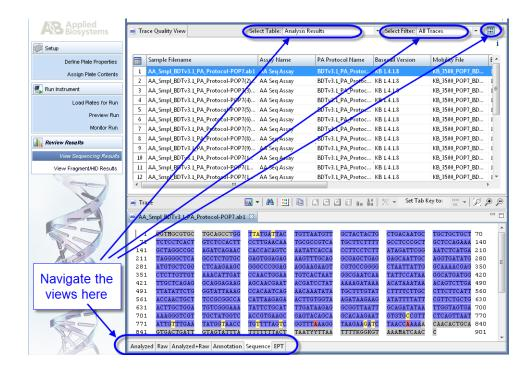
The 3500/3500xL analyzer display a progress indicator while it checks the level of consumables on the instrument.



**Monitor the run** Monitor the run by checking the status icons in the Injection Details section (Monitor Run screen).

Plate Name:	m Move Up in I	List 😻 Move Down in List 👔	🕽 Dølete   🕎 Duplicate 🚦	🖥 Re-Inject   🔟 Pause	Run 🔯 Resume Run 📗	🛛 Abort Injection 🛛 🔯 Terminate Injection I
AB Applied Biosystems		ntus: Connected Ime: Run 2009-04-39-11-52-26-9		: Administrator	Last Log Estimated Time Rei	in Time: 28-Apr-2009 12:43:51 PM naining: 03:05:54
Define Plate Properties Assign Plate Contents	<ul> <li>Injection List</li> <li>4 injections cre</li> </ul>	et Details eated = 4 in Plate A - 0 in Plate B				
💺 Run instrument	Injection	Type Assay	Instrument Protocol	Plate		'late B • • • • • • • • • • • • • • • • • •
Load Plates for Run	1 🔽	AA Assay 4 SeqScape	RapidSeq50_POP7xl_1	AA Seq Plate 4 Se		
Preview Run	2 🔼	AA Assay 4 SeqScape	RapidSeq50_POP7xl_1	AA Seq Plate 4 Se		2 3 4 5 6 7 8 9 10 11 12 🔺
	4	AA Assay 4 SeqScape AA Assay 4 SeqScape	RapidSeq50_POP7xl_1 RapidSeq50_POP7xl_1	AA Seq Plate 4 Se AA Seq Plate 4 Se		AA AA AA AA
Monitor Run					D AA	AA
2		/				AA
		/			F AA	AA
Review Results	$\cup$	/				

View sequencing results You can view the Sequencing Results in the 3500 Series Data Collection Software by going to the View Sequencing Results screen and selecting the tab of interest.



Confirm run<br/>completionWhen the run successfully transfers for downstream analysis, the AutoanalysisManager displays the project as successfully processed.

Edit Help				
neral SeqScape 2.7				
Job	Project	Arrival Date	Status	6
njection 2009-04-30-09-51-46-950 - Auto-Analysis_RG-BDTv3.1_PA_Protocol-POP7 - 2009-Apr-30-09-55-53	AA-Tester1-Project1	Apr 30, 2009 9:55:53 AM	Complete	Project successfully proces
< [	17			

You can now launch SeqScape® and review the analyzed project.

**Note:** For guidelines on reviewing data and results, see the *SeqScape® Software v2.7 Workflow Quick Reference Guide* (PN 4401740) or the *SeqScape® Software User Guide* (PN 4359442).

# Auto-analysis with MicroSeq<sup>®</sup> ID

For instructions detailing how to set up a MicroSeq<sup>®</sup> ID analysis protocol, see "Create a new MicroSeq<sup>®</sup> ID analysis protocol" on page 191. For installation information on setting up the MicroSeq<sup>®</sup> ID Software to work with the 3500 Series Data Collection Software, see the *MicroSeq<sup>®</sup> ID v2.2 Getting Started Guide*.

### Perform secondary analysis on fragment experiments

The Applied Biosystems 3500/3500xL Genetic Analyzers and 3500 Series Data Collection Software provide integration between the instrument and secondary fragment analysis software applications — specifically GeneMapper<sup>®</sup> Software v4.1 and GeneMapper<sup>®</sup> *ID-X* Software v1.1. Using auto-analysis, samples are loaded, fragment data is generated, and allele calling is performed according to the protocols assigned to the plates prior to the run.

Software	Purpose
GeneMapper <sup>®</sup>	A high-performing and versatile software package for all fragment analysis and genotyping applications.
GeneMapper <sup>®</sup> <i>ID-X</i>	A software for use in Human Identification testing (databasing, casework, and paternity applications) and used in conjunction with AmpF/STR kit and the 3500/3500xL analyzer.

### Auto-analyze projects in the fragment analysis software

Auto-analysis can only be performed on the same computer that collects the sample files, therefore GeneMapper<sup>®</sup> or GeneMapper<sup>®</sup> *ID-X* Software must be installed and configured with the 3500/3500xL analyzer on a Windows Vista<sup>®</sup> operating system. Secondary analysis occurs within the GeneMapper<sup>®</sup> or GeneMapper<sup>®</sup> *ID-X* Software.

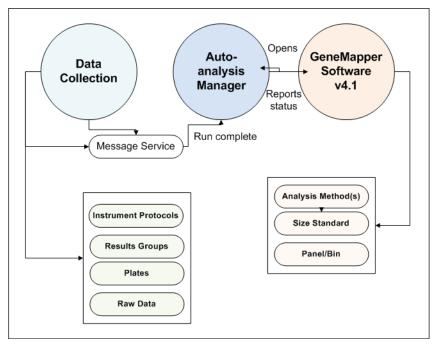
This procedure initially describes how to set up panels and bin sets in GeneMapper<sup>®</sup> Software v4.1 and then describes how to auto-analyze samples using the 3500 Series Data Collection Software. Once a run is complete, your data is seamlessly transferred into GeneMapper<sup>®</sup> for analyzing, processing and reporting.

**Note:** For detailed information on setting up a GeneMapper<sup>®</sup> *ID-X* analysis to autoanalyze in the 3500 Series Data Collection Software, see *GeneMapper<sup>®</sup> ID-X* v 1.1 *User Guide*.

# Set up an auto-analysis project in GeneMapper®

**IMPORTANT!** When using GeneMapper<sup>®</sup> Software to auto-analyze results data from the 3500/3500xL analyzer, you must have v4.1 installed on the *same* computer as the 3500 Series Data Collection Software.

The fragment analysis workflow for auto-analysis is summarized in this flow chart.



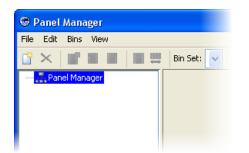
Set up a project in the secondary analysis software before starting a run on the 3500/3500xL analyzer. All analysis in GeneMapper<sup>®</sup> occurs in a project.

1. Open GeneMapper<sup>®</sup> v4.1 by double-clicking



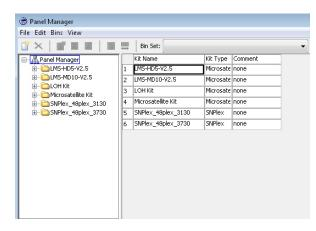
Specify a kit, a panel, and a bin set for the project

- 2. Click to open the Panel Manager.
- 3. Select the Panel Manager node (in the Navigation pane) to highlight.



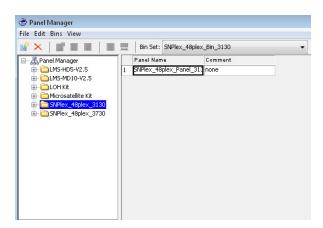
4. Select the panel, then **Import** to import a previously created kit folder with panel marker information.

Note: You have to import panels one by one; repeat this step for each panel.



5. Import the bin sets that are associated with the panels you just imported above. Click **Import** for each bin set.

Note: You have to associate a bin set to every panel that you imported.



6. Click **OK** to save and close the Panel Manager.

**Note:** For more information on how to create panels and bin sets, see the *GeneMapper® v4.1 Quick Reference Guide (PN 4362816)* or refer to the specific Getting Started Guide for your application.

# Create a new project

1. Click [] (GeneMapper Manager) to open the GeneMapper Manager.

2. Select the Analysis Method tab, then click New.

Name	Last Saved	Cwner	Instrument	Analysis Type	Description
SNPlex_Rules_3730	2004-09-21 11:39:02.0	gm	3730	SNPlex Rules 3730	Default SNPlex settings For 3730 or 3730xd
SNPlex_Rules_3130	2008-04-17 14:59:07.0	gm	3130	SNPlex Rules 3130	Default SNPlex settings for 3130 analysis
SNPlex_Model_3730	2004-09-21 11:38:48.0	gm	3730	SNPlex Model 3730	Default SNPlex settings for 3730 or 3730x1
SNaPshot Default	2003-07-25 18:10:08.0	gm		SNaPshot®	Factory Provided
Microsatellite Default	2005-06-08 13:50:00.0	gm		Microsatellite	Factory Provided
AFLP Default	2004-07-16 16:21:19.0	gm		AFLP	Factory Provided
<					D

3. Select the Analysis Method Type you want, then click **OK**.

New Analysis Method			<b>—</b>
Select analysis type:			
Microsatellite	🔘 OLA Analysis	AFLP	
SNaPshot®	🔘 SNPlex Model 3730		
SNPlex Rules 3730	SNPlex Rules 3130		
	OK Cancel		

4. Name your Analysis Method (General tab).

	or - Microsatellite eak Detector   Peak Quality   Quality Flags	
Analysis Method [	Description	
Name:	Test	
Description:		-
Instrument:	3500×L	
Analysis Type:	Microsatellite	

5. Select your Bin Set (Allele tab).

A

An	Analysis Method Editor - Microsatellite									
	General Allele Peak	Detector	Peak Quality	y Quality Fla	ags					
	Bin Set: None Marker Rept <mark>SWPlex_4</mark> Use rSNPlex_4 Values for dinucleo		_3730		e alle	•				
	values for ullitudeo	uue repea								
			Mono	Tri	Tetra	Penta	Hexa			
	Cut-off value		0.25	0.2	0.25	0.25	0.25			
	PlusA ratio		0.0	0.95	0.95	0.95	0.95			
	PlusA distance		0.0	1.6	1.6	1.6	1.6			
	Stutter ratio		0.0	0.95	0.15	0.15	0.15			
	Stutter distance	From To	0.0	0.0 3.5	0.0 4.5	0.0 5.5	0.0 6.5			

6. Select your Peak Detector Algorithm as Basic, Advanced or Classic (Peak Detector tab).

**IMPORTANT!** If you want to enable size standard normalization, you must select **Advanced**.

General Allele	Peak Detec	tor Peak	Quality	Quality Flag	js
Peak Detection #		Basic		-	
-Minimum Peak H	siant L	Basic Advanced			
		Elassic			
<ol> <li>Auto</li> </ol>	omatic				
🔘 Use	r specified (	rfu):			
	Green	Yellow	Red	Purple	Orange
Blue	0.000			50	50
Blue					1

7. Customize your Peak Quality and/or Quality Flag settings in the appropriate tab, then close the Analysis Method Editor.

Quality weights are b Quality Flag Settings	etween O	and 1.			
Quality Hug Socialitys					
Spectral Pull-up (SPL	0	0.5	Con	trol Concordance (CC	) 0.5
Broad Peak (BD)		0.5	Low	Peak Height (LPH)	0.5
Single Peak Artifact (SPA)		0.5	Off-	scale (OS)	0.5
Sharp Peak (SHP)		0.5	Pea	k Height Ratio (PHR)	0.5
Cross Talk (XTLK)		0.5	One Basepair Allele (OBA)		) 0.5
Out of Bin Allele (BIN)		0.8	Split Peak (SP)		0.5
PQV Thresholds					
	Pas	s Range:		Low Quality R	ange:
Sizing Quality:	From	0.75	to 1.0	From 0.0 to	D.25
Genotype Quality:	From	0.75	to 1.0	From 0.0 to	0.25

8. Click **OK** to save, then click **Done** to close GeneMapper<sup>®</sup>.

**IMPORTANT!** Close GeneMapper<sup>®</sup> v4.1 before performing the auto-analysis run on the 3500/3500xL analyzer.

# Set up a GeneMapper plate in the 3500 Series Data Collection Software

Set up a fragment analysis run in the 3500 Series Data Collection Software by assigning an Assay, a File Name Convention and a Results Group.

Start the 3500 Series Data Collection Software

- 1. Start the Auto-Analysis Manager before starting the 3500 Series Data Collection Software.
- 2. Start the 3500 Series Data Collection Software, then go Dashboard 🕨



3. Name your new plate.

late Details	
* Name:	Test Fragment Plate
* Number of Wells:	🖲 96 🔘 96-FastTube 🔘 384
* Plate Type:	Fragment 🔹
* Capillary Length:	50 🔻 cm
* Polymer:	POP7 -

- 4. Select the Number of Wells, Plate Type as Sequencing, Capillary Length and Polymer associated with this plate for the current run.
- 1. Check **Perform Auto-Analysis** (right side of the Plate Details section), to expand the Secondary Analysis section.

▼ Secondary Analysis		Perform Auto-Analysis 🕝
* Software Type:	GeneMapper 🔹	
* Software Location:	GeneMapper-3500-FVTEST 🔹	

2. Confirm GeneMapper auto-populates as the Software Type.

**Note:** If GeneMapper does *not* appear in the drop-down list under Software Type, check your installation. Secondary analysis software must be installed correctly before GeneMapper is automatically listed as a selection.

3. Confirm Your computer name auto-populates as the Software Location.

**IMPORTANT!** For auto-analysis to be successful, the secondary analysis protocol must match the software location set here.

- 4. Enter your Username and Password for auto-analysis access to the secondary analysis software.
- 5. Click Save Plate > Save to save your plate with these settings, then Assign Plate Contents to advance to the next screen.

analysis for secondary analysis

Specify auto-

**Assign plate contents** When assigning plate contents, you are assigning assays, file name conventions and results group to be associated with your auto-analysis.

Set up an assay 1. In the Assign Plate Contents screen, go to the Assays box and select either Create New Assay or Add From Library.

Assays		
	Actions	•
Add from Library		
Create New Assay		

2. Name your new assay in the Setup an Assay dialog box.

Note: (Optional) Select a color for this assay to display with in the Plate View.

3. Select an Instrument Protocol to apply to the assay.

**Note:** For more instruction on setting up an instrument protocol, see "Create a new instrument protocol" on page 165.

4. Select a Sizecalling Protocol to apply to the assay.

**Note:** For more instruction on setting up an instrument protocol, see "Create a new sizecalling protocol" on page 179.

5. Create a new fragment analysis protocol (GeneMapper Protocol) to apply to the samples by clicking **Create New**.

📜 Test Fragment F	Plate - Create New Ass	ay	X
Setup an Assa	у		
			0
* Assay Name:	Your Fragment Assa	y Docked	Color: Dark Blue 🔻
Application Type: Protocols Do you wish to		ment protocols to this assay? 💿 No 🛛 🔘 Yes	
* Instrument Pr	otocol:	FragmentAnalysis50_POP7xl_1	▼ Edit Create New
* Sizecalling Pr	otocol:	Fragment_Analysis_PA_Protocol	▼ Edit Create New
GeneMapper Pr	otocol:	Microsatellite SA Protocol SNPIex 3730 SA Protocol	Edit Create New
Close			bly to Plate Save to Library

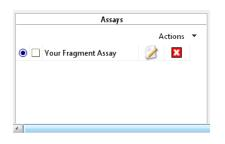
6. Name your new fragment analysis protocol and (optionally) enter a description.

7. Select the panel(s) you previously created in GeneMapper<sup>®</sup>, then click Apply to Assay.

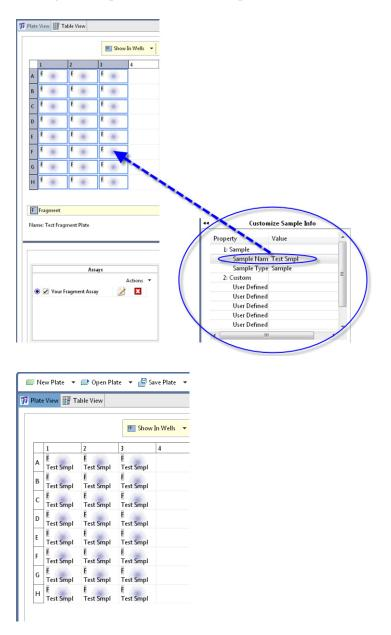
		condary Analysis Protocol		-
Setup a GeneM	lapper● Protocol			
				Ø
* Protocol Name:	Your Secondary Ana	alysis Protocol	C Locked	
Description:				
	Application Type:	Fragment Analysis	-	
* Seconda	ary Analysis Software:	GeneMapper	•	
* Secondary Analy	sis Software Instance:	GeneMapper-3500-FVTEST	•	
Properties				
	* Analysis Method:	Test	•	
	* Size Standard:	GS500(-35,-250,-340)ROX	•	
	Panel:	SNPlex_48plex_Panel_3130	•	
		SNPlex_48plex_Panel_3130	^	
		Panel11-HD5-V2.5 Panel02-HD5-V2.5		
				Apply to Assay Save to Library

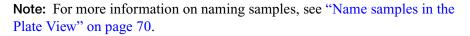
**Note:** For more instruction on setting up a secondary analysis protocol, see "Create a new fragment analysis protocol" on page 193.

- 8. Click **Close** when you are finished applying all the panels to the assay.
- 9. Click Apply to Plate, then close the Setup an Assay dialog box.



10. Name your samples by highlighting the number of wells in your plate and naming the sample in Customize Sample Info box.





Specify FNC and RG 1. Specify a File Name Convention (FNC) and a Results Group (RG) to associate with your project.

**Note:** You can create a FNC with the specimen name as a part of your sample file name.

- 2. Highlight the wells of your plate configuration (Plate View) and check the box next to the appropriate FNC to apply it to your project.
- 3. Repeat for the Results Group.

**Note:** For more information on setting up a FNC see "Create a new file name convention" on page 151. For more information on setting up a RG, see "Create a new results group" on page 156.

- 4. Click Save Plate > Save.
- 5. Click Link Plate for Run
- 6. Click Create Injection List, then click **OK** after the instrument performs its validations.

Start the autoanalysis run Click Start Run to begin your auto-analysis.

The 3500/3500xL analyzer displays a progress indicator while it checks the level of consumables on the instrument.

Progress Information	
Preparing Instrument Run	
Starting Instrument Run	
	Cancel

Confirm runWhen the run successfully transfers for downstream analysis, the AutoanalysiscompletionManager displays the project as successfully processed.

	Help GeneMapper v4.1 SeqScape 2.7						
	lapper Job Queue						
,	Job/Project	User	# of Samples	Arrival Date	Completed Date	Status	
	Injection 2009-05-04-17-34-59-945 - AutoAnalysisResultsGroup_Frag	gm	8	May 4, 2009 5:39:22 PM	May 4, 2009 5:47:06 PM	Complete	Analysis completed successfully
	Injection 2009-05-04-17-34-59-955 - AutoAnalysisResultsGroup_Frag	gm	8	May 4, 2009 5:40:41 PM	May 4, 2009 5:51:17 PM	Complete	Analysis completed successfully

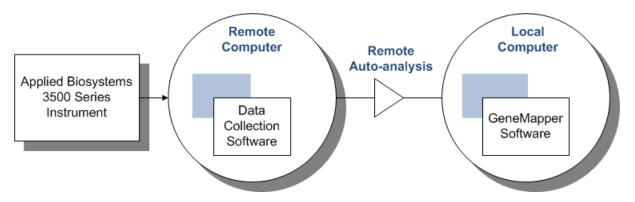
You can now launch GeneMapper<sup>®</sup> and review your analysis.

**Note:** For guidelines on reviewing fragment data and results, see the *GeneMapper*<sup>®</sup> v4.1 Quick Reference Guide (PN 4362816) or refer to the specific Getting Started Guide for your application.

### Auto-Analysis with GeneMapper® ID-X

For instructions detailing how to set up a GeneMapper<sup>®</sup> *ID-X* analysis protocol, see "Create a new HID analysis protocol" on page 195. For installation information on setting up the GeneMapper<sup>®</sup> *ID-X* Software v1.1 to work with the 3500 Series Data Collection Software, see the *GeneMapper<sup>®</sup> ID-X Software v1.1 User Guide*.

# Remote auto-analysis configuration



For remote auto-analysis, the 3500 Series Data Collection Software resides on the instrument computer and the GeneMapper<sup>®</sup> Software resides on a *different* computer.

In this configuration, you can set up both softwares so that GeneMapper<sup>®</sup>:

- connects to a remote computer running the 3500 Series Data Collection Software
- obtains sample files from the remote 3500 Series Data Collection Software database
- performs analysis of the generated sample files automatically

# Remote auto-analysis installation

Install the remote auto-analysis configuration when you want to auto-analyze data and you plan to connect to a separate computer running the 3500 Series Data Collection Software.

Installing GeneMapper<sup>®</sup> Software as a remote auto-analysis configuration requires that you:

- 1. Start the Data Collection services on the remote Data Collection computer.
- 2. Install GeneMapper<sup>®</sup> Software v4.1 on the local computer.

**IMPORTANT!** Before installing GeneMapper<sup>®</sup>, start the Data Collection services on the remote Data Collection computer.

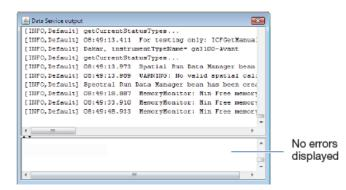
#### Start the 3500 data collection services

 On the Data Collection computer, select Start > All Programs > Applied Biosystems > Data Collection > Run Data Collection version 1.0.

Note: If the services do not start automatically, click Start All.

STOPPING POINT. Wait until all services have changed to green before continuing.

- 2. If the 3500 Series Data Collection Software requires a password, type the login name and password, then click **OK**.
- 3. Verify that Data Service started without errors:
  - a. In the Service Console, right-click the graphic next to each service listed and select **Show Console** to display the Data Service output window.
  - b. Verify that no errors are displayed, then close the Data Service dialog box.



- 4. Obtain the host name (full computer name):
  - a. Right-click **Computer** on the desktop, then select **Properties**.
  - b. Locate the full computer name. (You will need to enter the name when you install the GeneMapper<sup>®</sup> Software).

#### Windows Vista

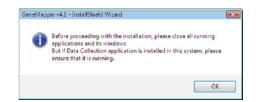
Computer name, domain, a	nd workgroup	settings
Computer name:	ServerName	_
Full computer name:	ServerName	)
Computer description:		-

c. Close the dialog box.

Install GeneMapper<sup>®</sup> v4.1 for remote auto-analysis 1. Insert the *GeneMapper*<sup>®</sup> v4.1 Software Full Installation DVD into the DVD drive to start the installer.

If the installer does not start automatically:

- a. Right-click Computer, then select Explore.
- b. Expand the DVD drive, then select the GeneMapper<sup>®</sup> v4.1 folder to display its contents.
- c. Double-click to start the installer.
- 2. Close all other applications and windows, then click **OK** to close this message:



- 3. In the Welcome window, click Next.
- 4. Review the installation requirements status, then click Next.
- 5. Select Remote Analysis for type of installation, then click Next.

GeneMapper	
GeneMapper® Software Version 4.1	Setup Type Phoose the setup type that best suits your needs.
A posted as	Remote auto-analysis option will instal the auto-analysis manager utility to automatically analyze data collected from Data Collection system. If you don't have a Data Eolection system available, you should choose stand-done option. Stand-alone Femote Autoanalysis
	< Back Next > Cancel

- 6. In the GeneMapper Client setup window, type the server name (full computer name) for the Data Collection computer (see step 4 on page 292), select ABI 3500, then click **Next**.
- 7. Read the release notes, then click Next.

**Note:** For other installation and configuration setup instructions, see Chapter 3 of the *GeneMapper*<sup>®</sup> v4.1 Installation and Administration Guide.

### Create a shared folder

Create a shared folder (Windows Vista<sup>®</sup>)

- 1. Select **Start** > **Computer**, then double-click the drive on which you want the shared folder to reside.
- 2. Select File ▶ New ▶ Folder.
- 3. Name the folder (for example: Remote\_Autoanalysis).
- 4. Right-click the new shared folder, then select Properties.
- 5. Select the Sharing tab, click Share.

📜 Remote_Aut	oanalysis Pro	operties		<b>.</b>	
General Shari	ng Security	Previous Versions	Customize		
Network File and Folder Sharing					
	emote_Autoan et Shared	alysis			
Network Path: Not Shared					
Share					

6. In the Choose people to share with dialog box, click the drop-down and select **Everyone (All users in this list)**.

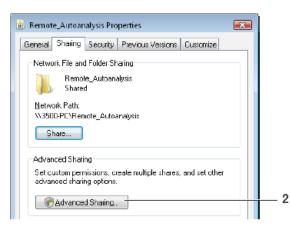
#### Choose people to share with

			- (
	-	Add	
3500-ADMIN			
3500-USER	) L	evel	
ABService			
Administrator			
veryone (All users in this list)			
reate a new user			
ell me about different ways to share in Windows.			
en me avoue amerene mejo eo anare ar runa onzi			

- 7. Click Add.
- 8. In the Permission Level column, change the value from Reader to Co-owner.
- 9. Click Share, then click Done and Close.

📔 Remote Aut	oanalysis (\\3500-P	0		
	emote_Autoanalys			

- 10. Click **OK**.
- Set security preferences for the shared folder
- 1. Right-click the shared folder, then select Properties.
- 2. Select the Sharing tab, then click Advanced Sharing.



3. On Advanced Sharing window, click Permissions.

Advanced Sharing	
Share this folder	
Settings	
Share name:	
Remote_Autoanclysis -	
Umt the number of smultaneous users to: 10 🚔	
Comments:	
	- 3
Permissions Caching	
OK Cancel Apply	

4. In the Permissions for the *<shared folder name>* dialog box, select the checkbox for **Full Control** (in the **Allow** column).

nare Permissions			
group or user names:			
🐍 3500-ADMIN (3500-PC\35	CO-AD MIN)		
	Add	Remove	
ermissions for 3500-ADMIN	Alow	Dery	
Full Control	<b>V</b>	1	
Change	$\checkmark$		
Read			

- 5. Click **OK** twice.
- 6. Click Close.

Turn off password protected sharing	<b>IMPORTANT!</b> Before starting Remote Auto-analysis, you must make sure that the password protected sharing settings on the Data Collection computer are turned off.
	<ol> <li>On the Data Collection computer, select Start &gt; Control Panel &gt; Network and Sharing Center.</li> </ol>

- 2. Click the expand button ( $\bigcirc$ ) for Password protected sharing,
- 3. Select Turn off password protected sharing.

Network and Sharing Co	enter	
3500-PC (This compute	applera.net	View full map
💐 applera.net (Public netwo	rk]	Customize
Access	Local and Internet	
Connection	AB Instrument - Motherboard NIC	View status
Sharing and Discovery Network discovery	© On	$\odot$
File sharing	On On	<ul> <li>Image: Second sec</li></ul>
Public folder sharing	On (password required)	$\odot$
Printer sharing	e Off	
Password protected sharing	© On	
	ected sharing	
3	5 OI	4

4. Click Apply.

# Set up the 3500 Series Data Collection Software v1.0

Complete autoanalysis setup on the Data Collection computer To complete your remote auto-analysis setup, you must create a new Results Group.

- 1. Select the Results Group node in the navigation pane.
- 2. Click New to open the Results Group editor.
- 3. Complete the selections in the General tab by:
  - a. Enter the new Results Group name.
  - b. Enter the Results Group owner.
  - c. (Optional) Enter the Results Group comment.
  - d. Check Results Group Entry Completed.
- 4. Complete the selections in the Analysis tab by:
  - a. Select the GeneMapper instance (GeneMapper + computer name) from the drop-down list.

b. Check **Do Autoanalysis**.

**Note:** If you plan to perform an auto-analysis for every Results Group Complete instead of each run individually, check **Results Group Entry Completed**.

- c. Enter the GeneMapper Login ID and password.
- 5. Complete the selections in the Destination tab by:
  - a. Check Use Custom Location.
  - b. Enter the Destination using the format:

\\*Remote analysis computer name*\*Shared folder name*, for example:

#### \\myPC\Remote\_Autoanalysis

- c. (For remote auto-analysis specifically) Establish a connection with the remote analysis computer by:
  - Select Start > Run.
  - Enter the destination path, then click **OK**.
  - Click **Test** to test the Location path name connection.

**STOPPING POINT.** If the test **Passes**, the message displays "Path Name test successful." If the test **Fails**, the message displays "Could not make the connection. Please check that the Path Name is correct." In this case, click **Browse**, then select the correct location.

If you encounter any unforeseen and potentially hazardous event while operating the instrument, turn off the power switch, unplug the instrument, and call your Applied Biosystems representative.

**IMPORTANT!** See the Safety appendix for instrumentation and chemical safety information and guidelines.

### Instrument troubleshooting

Symptom	Possible cause	Action	
Amber light (blinking)	Run paused	Resume run	
	Door open	Close the instrument door	
	Run failure that doesn't require restart of instrument	Conduct another run	
Instrument status light is blinking red	Instrument error	1. Power off the instrument.	
		2. Power on the instrument.	
		3. Restart the computer.	
"An error has been detected from the instrument."	Instrument monitor circuit failure	Restart the computer	
<complex-block></complex-block>	One or more of the services are stopped.	Right-click the status icon, then select Services. If any item does not display a checkmark, click the item to start the service. Services	
"Unable to transmit measurement data. Internal data buffer overflow."	Communications error.	Restart instrument and computer.	
Electric discharge message during	The ABC buffer may be low.	Replace the ABC.	
runs.		Ensure that the ABC is being replaced per 3500 Series Data Collection Software notifications.	

# Spatial calibration troubleshooting

Symptom	Possible cause	Action
"Start" Spatial Calibration button is disabled.	Communication failure between the Data Collection Software and instrument	Restart instrument and computer. Check the NIC cable connection.
Unusual peaks or a flat line for the spatial calibration.	Improper installation of the detection cell: Detection cell on the array is not properly seated.	<ul> <li>Unistall, then re-install the array: Reinstall the detection cell to reposition and make sure it fits in the proper position. If the calibration fails again:</li> <li>1. Fill the capillaries with polymer.</li> <li>2. Repeat the spatial calibration.</li> </ul>
	The instrument may need more time to reach stability. An unstable instrument can cause a flat line with no peaks in the spatial view.	Repeat the spatial calibration.
	Broken capillary resulting in a bad array fill.	Check for a broken capillary, particularly in the detection cell area. If necessary, replace the capillary array using the Wizard.
Persistently bad spatial calibration results.	Bad capillary array.	Replace the capillary array, and then repeat the calibration. Call your Applied Biosystems representative if the results do not improve.
"Spatial Calibration Error" message. The instrument cannot perform Spatial Calibration with Array fill.	Conditioning reagent is installed.	Replace the Conditioning reagent with an appropriate Polymer.

# Spectral calibration troubleshooting

Symptom	Possible cause	Action
No signal	Incorrect preparation of sample	Replace samples with fresh samples prepared with fresh Hi-Di <sup>™</sup> Formamide
	Bubbles in sample wells	Centrifuge samples to remove bubbles
	The capillary tips may not be touching the samples.	Check the volume of your samples. If no results, call your Applied Biosystems representative.
	The capillary tips may be hitting the bottom of the wells. Autosampler not correctly aligned.	Call your Applied Biosystems representative.
If the spectral calibration fails, or if a message displays "No spectral files found."	Blocked capillary	Refill the capillary array. You may have to install a fresh array or consider that capillary non-usable for purposes of planning your runs.
	Incorrect chemistry file, dye set, and/or run module selected.	Correct the files and rerun the calibration.
	Insufficient filling of array.	Check for broken capillaries and refill the capillary array.
	Expired matrix standards or old reagents.	Check the expiration date and storage conditions of the matrix standards and/or reagents. If necessary, replace with a fresh lot.
Data Error - One or more peaks fall below the minimum required amplitude of 750.	One or more peaks fall below the minimum required amplitude of 750.	Rerun the spectral standards, and if necessary, increase the amount of spectral standard added.
Spikes in the data or "Bad dye order detected" error message.	Expired polymer.	Replace the polymer with a fresh lot using the Replenish Polymer Wizard.
	Bubbles in the polymer system.	Select the Bubble Remove Wizard to clear the bubbles.
	Possible contaminant or crystal deposits in the polymer.	Properly bring the polymer to room temperature; do not heat. Replace the polymer if it has expired.
Elevated baseline.	Poor spectral calibration.	Perform new spectral calibration.
Spectral calibration history does not display previously run calibration.	If you change polymer type, spectral calibrations for the original polymer type are not retained.	No action.
Pull-down (mirror image) peaks	The first time you perform a spectral calibration (for each dye set) after installing a new capillary array, you may notice pull-down peaks (or mirror image peaks). These pull-down peaks will eventually correct themselves once the run completes.	No action.

# Sequencing install standard troubleshooting

Symptom	Possible cause	Action
No signal	Incorrect preparation of sample	Replace samples with fresh samples prepared with fresh Hi-Di <sup>™</sup> Formamide.
	Bubbles in sample wells	Centrifuge samples to remove bubbles.
	The capillary tips may not be touching the samples.	Check the volume of your samples. If no results, call your Applied Biosystems representative.
	The capillary tips may be hitting the bottom of the wells. Autosampler not correctly aligned.	Call your Applied Biosystems representative.
If the Sequencing install standard (Performance check) fails. Fail capillary • If more than one failed capillary (for 8-capillary). • If more than three failed capillary (for 24-capillary). Accept button is not active, but Reject button is active.	Blocked capillary	Refill the capillary array. You may have to install a fresh array or consider that capillary non-usable for purposes of planning your runs.
	Incorrect chemistry file, dye set, and/or run module selected.	Correct the files and rerun the calibration.
	Insufficient filling of array.	Check for broken capillaries and refill the capillary array.
	Expired matrix standards or old reagents.	Check the expiration date and storage conditions of the matrix standards and/or reagents. If necessary, replace with a fresh lot.
	Expired polymer.	Replace the polymer with a fresh lot using the Replenish Polymer Wizard.
	Bubbles in the polymer system.	Select the Bubble Remove Wizard to clear the bubbles.
	Possible contaminant or crystal deposits in the polymer.	Properly bring the polymer to room temperature; do not heat. Replace the polymer if it has expired.

# Fragment/HID install standard troubleshooting

Symptom	Possible cause	Action Select all dyes, then generate the report.				
Fragment/HID report contains blank pages or incomplete information.	All dyes are not selected before you generate the report.					
No signal	Incorrect preparation of sample	Replace samples with fresh samples prepared with fresh Hi-Di <sup>™</sup> Formamide				
	Bubbles in sample wells	Centrifuge samples to remove bubbles.				
	The capillary tips may not be touching the samples.	Check the volume of your samples. If no results, call your Applied Biosystems representative.				
	The capillary tips may be hitting the bottom of the wells. Autosampler not correctly aligned.	Call your Applied Biosystems representative.				
If the Fragment/HID install standard (Performance check) fails.	Blocked capillary	Refill capillary array. You may have to install a fresh array or consider that capillary non-usable for purposes of planning your runs.				
	Insufficient filling of array.	Check for broken capillaries and refill the capillary array.				
	Expired matrix standards or old reagents.	Check the expiration date and storage conditions of the matrix standards and/or reagents. If necessary, replace with a fresh lot.				
	Expired polymer.	Replace the polymer with a fresh lot using the Replenish Polymer Wizard.				
	Bubbles in the polymer system.	Select the Bubble Remove Wizard to clear the bubbles.				
	Possible contaminant or crystal deposits in the polymer.	Properly bring the polymer to room temperature; do not heat. Replace the polymer if it has expired.				

# Anode buffer container troubleshooting

#### Also see "Data/electropherogram troubleshooting" on page 306.

Symptom Possible cause Action					
Buffer below fill line (inadequate amount of buffer).	Ensure that buffer level is at or above the fill line. Do not use if buffer level is too low or seal has been compromised.				
	Buffer below fill line (inadequate				

# Cathode buffer container troubleshooting

Also see "Data/electropherogram trou	ibleshooting" on page 306.
--------------------------------------	----------------------------

Symptom	Possible cause	Action
Electrophoresis failure.	Buffer below fill line (inadequate amount of buffer).	Ensure that buffer level is at or above the fill line.
		Do not use if buffer level is too low or seal has been compromised.

# **RFID** troubleshooting

Symptom	Possible cause	Action
Unable to read RFID information. "Failure to Read from RFID tag"	Consumable package is improperly installed or defective label.	Ensure that the RFID label is not visibly damaged and consumable package is
	Polymer/Conditioning reagent	properly installed.
	pouch mis-oriented.	Ensure that label is close, and parallel, to the instrument.
		Reposition or re-install consumable, and click <b>Refresh</b> on the dashboard.
		If no results, restart the instrument and the computer.
		If no results, install a new consumable (if available), and call your Applied Biosystems representative for a replacement.

# Link a plate troubleshooting

Symptom	Possible cause	Action
Plate does not link.	Spatial/Spectral calibration was not performed	<ol> <li>Perform spatial calibration.</li> <li>Relink the plate(s).</li> </ol>
Plate was linked, but now it is unlinked.	If you access the Load Plates for Run screen from the navigation pane, a plate may not be linked (indicated by the active Link button).	Access the Load Plates for Run screen from the navigation pane and click <b>Link Plate</b> .
"No plate in position A" message.	You physically loaded plate in position B (plate B position) and try to link plate.	Click <b>Link Plates</b> and link the plate directly to position B (plate B position). Follow the prompts.
"No plate detected" message	The plate is in position B.	Place the plate in position A.

# How to search and use the log files

The 3500 Series Data Collection Software generates the following log files that you view using a text editor such as Wordpad:

- 3500UsageStatistics.txt–Provides a summary of the number of plates run, as well as number of run types (sequencing, fragment, and HID).
   Stored in: x:\Applied Biosystems\3500\UsageData
- **3500ConsumableUpdates.txt**–Provides a summary of consumables installation information and dates.

Stored in: D:\Applied Biosystems\3500\LogFiles

# View instrument sensor details

Click **View Instrument Sensor Details** in the Dashboard to display instrument information.

View Instrument Sensor Details

Run status of the instrument is displayed while a run is in progress.

Sensor States	Sensor Values											
Laser: On	EP Voltage (kV)		CEP Curre	nt (µA) ——		Laser	Powe	er (mW)—		Laser C	urrent (mA) —	
EP: On	19.7 🕨 💶 🕇 19.7	20.0		010	500.0		-	10	20.0		212	3000.0
Oven: Off				1 1			_	- Suites			1	
Oven Door: Open		15.0			375.0	13.4		iat:15.3	15.0			2250.0
Messages		10.0			250.0		-	-	10.0			1500.0
		5.0	119.8		125.0		-	-	5.0			750.0
	-   -	0.0		-   -	0.0		-	-	0.0	12.1	⊾l-	0.0
	Oven Temp (°C)		CCD Can	nera Temp (	°C)	Ambie	ent Te	emp (°C)-		Detectio	on Cell Temp (	PC)
	111	70.0		111	0.0		_	12	40.0		111	70.0
	53.5 ⊾ <b>4t55.</b> 4	60.0		1 1	-4.0		_	-	35.0			60.0
	53.5 E	50.0					_	-	30.0			50.0
		40.0	11.0	1 1	-8.0	20.9	-	-	25.0			40.0
		30.0	-11.2		-12.0	20.9		1	20.0		 ■ <b>4t:30.0</b>	
		20.0			-16.0		-	-	15.0 10.0	23.5		20.0
	-   -	10.0		-   -	-20.0		-	-	5.0		-   -	10.0

Figure 37 Instrument sensor details

# Data/electropherogram troubleshooting

Symptom	Possible cause	Action					
Signal too high.	Sample concentration is too high.	Dilute the sample.					
		Decrease the injection time.					
	Too much DNA added to the reaction, resulting in uneven signal distribution.	Optimize reaction conditions.					
No signal.	Failed reaction.	Repeat reaction.					
	Blocked capillary.	Refill capillary array. You may have to install a fresh array or consider that capillary non-usable for purposes of planning your runs.					
	Bent capillary array tips.	Replace the capillary array.					
	Cracked or broken capillary array.	Visually inspect the capillary array, including the detector window area for signs of breakage.					
Low signal strength.	Degraded Formamide.	Use a fresh aliquot of Hi-Di <sup>™</sup> Formamide.					
	Not enough sample: Pipetting error.	Increase the amount of DNA added.					
	Sample has high salt concentration.	Dilute with distilled or deionized water.					
		Desalt using a column purification method.					
	Insufficient mixing.	Vortex the sample thoroughly, and then centrifuge the tube to condense the sample to the bottom of the tube.					
	Weak amplification of DNA.	Reamplify the DNA.					
		Check DNA quality.					
	Autosampler out of calibration.	Check the volume of your samples. If still low signal strength, call your Applied Biosystems representative.					
Elevated baseline.	Possible contaminant in the polymer path.	Use the conditioning reagent for washing the polymer pump.					
	Possible contaminant or crystal deposits in the polymer.	Bring the polymer to room temperature.					
		Replace the polymer if it has expired.					
	Poor spectral calibration.	Perform new spectral calibration.					
Loss of resolution.	Too much sample injected.	Dilute the sample and re-inject.					
	Poor quality water.	Use distilled or deionized water.					
	Degraded polymer.	Use a fresh supply of polymer.					
	Capillary array used for more than 160 injections.	Replace with new capillary array.					
	Degraded formamide.	Prepare fresh Hi-Di <sup>™</sup> Formamide and re-prepare samples.					
	High salt concentration in samples.	Use a recommended protocol for salt removal.					
		Dilute salts with water.					

Symptom	Possible cause	Action
Poor resolution in some capillaries.	Insufficient filling of capillary array.	Refill the capillary array and look for polymer leakage. If problem persists, call your Applied Biosystems representative.
		Re-inject the same samples.
	Poor quality samples.	Check the sample preparation.
	Leak in system.	Tighten the connectors and array lever
No current.	Not enough buffer in ABC.	Ensure that the buffer is filled up to the fill line.
	Bubble(s) present in the lower polymer block and/or the array	Pause the run and inspect for bubbles hidden in the tubing connectors.
	and/or channels.	Select the Bubble Remove Wizard to remove the bubbles.
Elevated current.	Degraded polymer.	Open fresh supply of polymer and use Replenish Polymer Wizard.
	Arcing in the lower polymer block.	Inspect the lower polymer block for discoloration or damage.
		Replace the lower polymer block if necessary.
Fluctuating current.	Bubble in polymer block.	Pause run and inspect for bubbles hidden in the tubing connectors.
		Select Bubble Remove Wizard to remove the bubbles.
	A slow leak may be present in the	Check polymer blocks for leaks.
	system.	Tighten all fittings.
	Not enough buffer in ABC.	Ensure that the buffer is filled up to the fill line.
	Arcing	Check for moisture in and around the septa, the CBC, the oven, and the autosampler.
Poor performance of capillary array used for fewer than 100 runs.	Poor quality samples, possible cleanup problems.	Desalt samples using a recommended purification protocol.
	Poor quality formamide.	Prepare fresh Hi-Di <sup>™</sup> Formamide and re-prepare samples.
	Leak in system.	Tighten the connectors and array leve
Migration time becomes	Leak in system.	Tighten the connectors and array leve
progressively slower.	Improper filling of the system with polymer.	Polymer delivery pump may need to be serviced.
		If the issue persists, call your Applied Biosystems representative.
Migration time becomes progressively faster.	Water in polymer system, resulting in diluted polymer.	Use Bubble Remove Wizard to add polymer to system.
	Buffer valve leakage.	Check the Buffer-Pin Valve and see if i closes correctly.

Symptom	Possible cause	Action
Extra peaks in the electropherogram.	Data off scale.	Dilute the sample and re-inject the sample.
	Possible contaminant in sample.	Re-amplify the DNA.
	Sample re-naturation.	Heat-denature the sample in good quality formamide and immediately place on ice.
Peaks exhibit a shoulder in GeneMapper <sup>®</sup> <i>ID-X</i> Software applications.	Sample re-naturation.	Heat-denature the sample in good quality formamide and immediately place on ice.
Error messages:	Bubbles in the polymer system.	Select the Bubble Remove Wizard to clear bubbles.
polymer delivery"	Leak in the polymer system.	Check for evidence of leaks.
<ul> <li>"Leak detected during bubble compression"</li> <li>The run aborts.</li> </ul>		If polymer leak occurred, conduct a water wash and wash the pump trap using the cleaning kit supplied.
	Buffer valve leakage.	Check the Buffer-Pin Valve and see if closes correctly.
		Clean the Buffer-Pin Valve.
		Ensure that the maintenance schedul is followed per 3500 Series Data Collection Software notifications.
	Filling the array during install array.	Run Fill the Array with fresh Polymer wizard, or run Change Polymer Type wizard.
Detection cell stuck.	Improperly placed detection cell.	To loosen the detection cell:
It is difficult to remove when changing the capillary array.		<ol> <li>Undo the array lever and pull the polymer block towards you to first notch.</li> </ol>
		<ol> <li>Hold both sides of the capillary array around the detection cell are and apply gentle pressure equally on both sides.</li> </ol>
		3. Release.
Electrophoresis current is unstable	Bubbles in the polymer system.	Select the Bubble Remove Wizard to clear bubbles.
Electrophoresis failure.	Buffer below fill line.	Ensure that the buffer has not split in the overflow. If so, move the buffer back to main reservoir.

## Dashboard troubleshooting

Symptom	Possible Cause	Action
The Days Remaining value for buffer/polymer does not automatically update.	The Days Remaining for buffers updates only when you click <b>Refresh</b> or <b>Start A Run</b> .	As part of daily startup, click <b>Refresh</b> to update buffer status.

## Load plate troubleshooting

Symptom	Possible Cause	Action
Pre-run validation check does not lisplay a date for a consumable.	The software does not display a date if it is identical to the preceding date. In the example below, the installation and recommended replacement dates for cathode buffer are identical to the dates for anode buffer.	No action.
EInstrument Run  Pre-run validation check failed		
Reason -	OK << Details	
Reason - Consumable: Anode Buffer has exceeded recomme Installation date: Jan J, 2009 Recommended replacement: Jan 8, 2009 Consumable: Cathode Buffer has exceeded recomm	nded time on instrument	
Consumable: Anode Buffer has exceeded recomme Installation date: Jan 1, 2009 Recommended replacement: Jan 8, 2009	nded time on instrument	Click <b>Details</b> to determine the cause of the error.

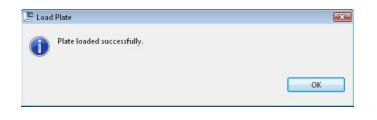
No sample name has been assigned to the following well: G01 H01 G02 H02 G03 44. To proceed, please assign a sample name to the wells Incompatible Polymer type Configured on plate: POP7 Currently installed: POP4 Incompatible Capillary length Configured on plate: 50 cm

OK

<< Det de

5

Symptom	Possible Cause	Action
"No plate in position A" message.	You physically loaded plate in position B (plate B position) and try to link plate.	Click <b>Link Plates</b> and link the plate directly to position B (plate B position).
		Follow the prompts.
"No plate detected" message	The plate is in position B.	Place the plate in position A
"Fragment performance check is	Running fragment modules after loading the plate.	Change polymer to POP-7 <sup>™</sup> .
required" message.		Run fragment analysis performance check.
"Sequencing performance check is	Running sequencing modules (POP-6 <sup>™</sup> ) after loading the plate.	Change polymer to POP-7 <sup>™</sup> .
required" message.		Run sequencing performance check.
After loading the plate.		
Load plate or Load Plate for Run message.	Performance issues.	Click <b>OK</b> and follow the prompts.



LUa	d Plate for Run
5	One or more positions on the instrument stacker are already occupied.
5)	Current injection list (if any) will be cleared.
	Please select one of the options below.
) Pla	ce this plate onto the stacker in position A, replacing the plate that is already in that position
Loa	nd this plate onto the stacker in position B
	OK Cancel

## Monitor run troubleshooting

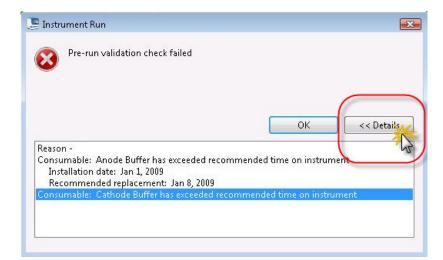
Symptom	Possible Cause	Action
Re-inject button is dimmed when you select an injection	Injection contains samples with assays that specify more than one instrument protocol.	Select in the injection list the injection with the instrument protocol of interest, select in the array view the capillary that corresponds to the well of interest, then click Re-inject.
The instrument run goes into pause state, unexpectedly.	RFID read/write process.	Check the Dashboard. Conduct an RFID refresh, if it does not refresh, restart both the computer and instrument.
Start run does not respond	The instrument has not initialized.	It takes, approximately, 10 seconds for the instrument to initialize after the instrument door is closed. Do not start a run until the instrument status light is green.
"Fragment performance check is required" message.	Running fragment modules after loading the plate.	Change polymer to POP-7 <sup>™</sup> . Run fragment analysis performance check.
"Sequencing performance check is required" message. After loading the plate.	Running sequencing modules (POP-6 <sup>™</sup> ) after loading the plate.	Change polymer to POP-7 <sup>™</sup> . Run sequencing performance check.

## Review results troubleshooting

Symptom	Possible Cause	Action
Sample files are not displayed when imported.	You imported (.hid) files and you did not click HID Samples.	Click HID Samples.
Peaks are not labeled when you access the screen.	Labels are not automatically applied.	See "Label peaks" on page 93.
x and y scaling plot settings are not applied when you click Apply.	Scaling settings are applied only when you click Zoom.	Click <b>Zoom</b> .
The sizing quality result reported in the 3500 Series Data Collection Software differs from the sizing	You imported (.fsa) files instead of (.hid) files into the GeneMapper <sup>®</sup> <i>ID-X</i> Software.	No action.
quality result for reported in the GeneMapper <sup>®</sup> <i>ID-X</i> Software.	The 3500 Series Data Collection Software <i>does not</i> consider the presence of broad peaks when determining sizing quality for fragment analysis data, therefore the sizing quality result reported in the 3500 Series Data Collection Software will differ from the sizing quality result reported in the GeneMapper <sup>®</sup> <i>ID-X</i> Software, which considers broad peaks in sizing quality.	

### Review error message details

Error messages in the 3500 Series Data Collection Software include a Details button. Click Details to display more information about an error message.



## Audit troubleshooting

Symptom	Possible Cause	Action
"Export did not complete successfully"	You exported records for samples that are not in their original location (samples have been deleted or moved).	Return sample data files to their original location, then export again.

## Electronic signature troubleshooting

Symptom	Possible Cause	Action
Electronic signature prompt is displayed when you edit sample comments.	Electronic signature prompt is displayed for sample comments, regardless of the electronic signature setting.	No action.

## Manual commands troubleshooting

Symptom	Possible Cause	Action
When you select Tools > Manual Commands, Set defined command for Consumables, then select a Read Command, the information displayed is not readable.	The feedback from Consumables Read Tag commands does not display valid information.	Refer to the Dashboard for consumables RFID tag information.

## Miscellaneous

Symptom	Possible Cause	Action
Polymer crystals on the Buffer-Pin	Buffer valve leakage.	Clean the Buffer-Pin Valve.
Valve.		Ensure that the maintenance schedule is followed per 3500 Series Data Collection Software notifications.
Fluid does not move through the pump and into the ABC from polymer or conditioning pouch.	Not applicable.	Call your Applied Biosystems representative.
Electric discharge message during	ABC may be low.	Replace the ABC.
runs.		Ensure that the ABC is being replaced per 3500 Series Data Collection Software notifications.
Leak detected during bubble compression during run or while filling the array.	Leak in system.	Run the Bubble Removal wizard. Ensure that there are no bubbles in the pump.
		If problem persists, use conditioning pouch for water wash. Use Replenish Polymer wizard to fill pump and array with polymer.
Only some injections, from a series of injections, are completed.	3500 Series Data Collection Software never moves on to the next injection.	Check connection between the instrument and computer and restart
"Injection failed" message.	Capillary RFID cannot be read.	both the instrument and computer.
After some of the injections complete.		Set up the injections again and started the runs.
When you click <b>Refresh</b> on the dashboard, and consumables information is listed as "Unknown."	Connection between the computer and instrument.	
"Instrument is not connected" message.	-	
After you start 3500 Series Data Collection Software.		
"Internal buffer data overflow" message.		

### Reset the instrument

Reset the instrument when:

- There is a fatal error as indicated by the red status light
- The instrument does not respond to the Data Collection software

Reset with the Reset button

1. Shut down the computer.

- 2. Close the instrument doors.
- 3. Reset the instrument with the Reset button, as shown.

**Note:** The Reset button is accessible through a small hole to the left of the Tray button.



#### Reset button

Reset by powering down

1. Shut down the computer.

- 2. Close the instrument doors.
- 3. Power off the instrument by pressing the on/off button on the front of the instrument.

4. Power on the instrument and wait until indicator light turns solid green.

- 5. Power on the computer.
- 6. Launch the Data Collection software (Service Console applications start automatically).

**IMPORTANT!** Wait until the computer has completely restarted before proceeding.

## Instrumentation safety

#### Symbols on instruments

Electrical symbols on instruments The following table describes the electrical symbols that may be displayed on Applied Biosystems instruments.

Symbol	Description	
-ÿ;-	Light switch.	
	Indicates the <b>On</b> position of the circuit breaker.	
0	Indicates the <b>Off</b> position of the circuit breaker.	
G	Indicates a standby switch by which the instrument is switched on to the <b>Standby</b> condition. Hazardous voltage may be present if this switch is on standby.	
0	Indicates the <b>On/Off</b> position of a push-push main power switch.	
Ŧ	Indicates a terminal that may be connected to the signal ground reference of another instrument. This is not a protected ground terminal.	
	Indicates a protective grounding terminal that must be connected to earth ground before any other electrical connections are made to the instrument.	
~	Indicates a terminal that can receive or supply alternating current or voltage.	

**Safety symbols** The following table describes the safety symbols that may be displayed on Applied Biosystems instruments. Each symbol may appear by itself or with text that explains the relevant hazard (see "Safety labels on instruments" on page 317). These safety symbols may also appear next to DANGERS, WARNINGS, and CAUTIONS that occur in the text of this and other product-support documents.

315

Symbol	Description	
$\triangle$	Indicates that you should consult the manual for further information and to proceed with appropriate caution.	
4	Indicates the presence of an electrical shock hazard and to proceed wit appropriate caution.	
	Indicates the presence of a hot surface or other high-temperature hazard and to proceed with appropriate caution.	
	Indicates the presence of a laser inside the instrument and to proceed with appropriate caution.	
	Indicates the presence of a biological hazard and to proceed with appropriate caution.	
	Indicates the presence of sharp object and piercing injury and to proceed with appropriate caution.	

#### Environmental symbols on instruments

The following symbol applies to all Applied Biosystems electrical and electronic products placed on the European market after August 13, 2005.

Symbol	Description
	<b>Do not dispose of this product as unsorted municipal waste.</b> Follow local municipal waste ordinances for proper disposal provisions to reduce the environmental impact of waste electrical and electronic equipment (WEEE).
	<b>European Union customers:</b> Call your local Applied Biosystems Customer Service office for equipmen pick-up and recycling. See <b>www.appliedbiosystems.com</b> for a list of customer service offices in the European Union.

### Safety labels on instruments

The Applied Biosystems 3500/3500xL Genetic Analyzers contain warnings at the locations shown below:

#### Locations of laser warnings

On the detection cell as shown below.



#### General instrument safety

WARNING! PHYSICAL INJURY HAZARD. Use this product only as specified in this document. Using this instrument in a manner not specified by Applied Biosystems may result in personal injury or damage to the instrument.
WARNING! PHYSICAL INJURY HAZARD. Using the instrument in a manner not specified by Applied Biosystems may result in personal injury or damage to the instrument.
<b>CAUTION! PHYSICAL INJURY HAZARD.</b> The instrument is to be moved and positioned only by the personnel or vendor specified in the applicable site preparation guide. If you decide to lift or move the instrument after it has been installed, do not attempt to lift or move the instrument without the assistance of others, the use of appropriate moving equipment, and proper lifting techniques. Improper lifting can cause painful and permanent back injury. Depending on the weight, moving or lifting an instrument may require two or more persons.
WARNING! Do not attempt to lift or move the computer or the monitor without the assistance of others. Depending on the weight of the computer and/or the monitor, moving them may require two or more people. Things to consider before lifting the computer and/or the monitor:

Things to consider before lifting the computer and/or the monitor:

 $\langle \cdot \rangle$ 

	• Make sure that you have a secure, comfortable grip on the computer or the monitor when lifting.
	• Make sure that the path from where the object is to where it is being moved is clear of obstructions.
	• Do not lift an object and twist your torso at the same time.
	• Keep your spine in a good neutral position while lifting with your legs.
	• Participants should coordinate lift and move intentions with each other before lifting and carrying.
	• Instead of lifting the object from the packing box, carefully tilt the box on its side and hold it stationary while someone slides the contents out of the box.
Operating the	Ensure that anyone who operates the instrument has:
instrument	• Received instructions in both general safety practices for laboratories and specific safety practices for the instrument.
	<ul> <li>Read and understood all applicable Material Safety Data Sheets (MSDSs). See "About MSDSs" on page 329.</li> </ul>
Cleaning or decontaminating	<b>CAUTION!</b> Before using a cleaning or decontamination method other than those recommended by the manufacturer, verify with the manufacturer that the

proposed method will not damage the equipment.

#### Physical hazard safety

the instrument

Moving parts

**WARNING! PHYSICAL INJURY HAZARD.** Keep hands clear of moving parts while operating the instrument. Disconnect power before servicing the instrument.

Solvents and pressurized fluids



**WARNING!** PHYSICAL INJURY HAZARD. Always wear eye protection when working with solvents or any pressurized fluids.

#### **Electrical safety**

WARNING! ELECTRICAL SHOCK HAZARD. Severe electrical shock ′5` can result from operating the 3500/3500xL analyzers without its instrument panels in place. Do not remove instrument panels. High-voltage contacts are exposed when instrument panels are removed from the instrument.

Power

WARNING! ELECTRICAL HAZARD. Grounding circuit continuity is required for the safe operation of equipment. Never operate equipment with the grounding conductor disconnected.

WARNING! ELECTRICAL HAZARD. Use properly configured and approved line cords for the voltage supply in your facility.



WARNING! ELECTRICAL HAZARD. Plug the system into a properly grounded receptacle with adequate current capacity.

Overvoltage The 3500/3500xL analyzers system has an installation (overvoltage) category of II, rating and is classified as portable equipment.

#### Laser safety

The 3500 or 3500xL analyzer uses a a solid state laser. The laser specifications are: Laser classification

- Wavelength 505nm ٠
  - Output power 20mW ٠

The LED specifications are:

- Emitting color Natural White ٠
- Luminous Intensity 250 Cd ٠

Under normal operating conditions, the instrument is categorized as a Class I laser product. When safety interlocks are disabled during certain servicing procedures, the laser can cause permanent eye damage, and, therefore, is classified under those conditions as a Class 3B laser.



CAUTION! LASER. Use of controls or adjustments or performance of procedures other than those specified herein may result in hazardous radiation exposure.

Laser safety requirements	<ul> <li>To ensure safe laser operation:</li> <li>The system must be installed and maintained by an Applied Biosystems Technical Representative.</li> <li>All instrument panels must be in place on the instrument while the instrument is operating. When all panels are installed, there is no detectable radiation present and the instrument is Class I. If any panel is removed when the laser is operating (during service with safety interlocks disabled), you may be exposed to laser emissions in excess of the Class 3B rating.</li> </ul>
	• Do not remove safety labels or disable safety interlocks.
Additional laser safety	Refer to the user documentation provided with the laser for additional information on government and industry safety regulations.
information	Also, note the laser warnings provided in "Safety labels on instruments" on page 317.
	WARNING! LASER HAZARD. Lasers can burn the retina, causing permanent blind spots. Never look directly into the laser beam. Remove jewelry and other items that can reflect the beam into your eyes. Do not remove the instrument top or front panels. Wear proper eye protection and post a laser warning sign at the entrance to the laboratory if the top or front panels

are removed for service.

#### Bar code scanner laser safety

Using a bar code scanner is optional.

Laser classification

Laser safety requirements

Class 2 (II) lasers are low-power, visible-light lasers that can damage the eyes. Never look directly into the laser beam. The scanner is designed to prevent human access to harmful levels of laser light during normal operation, user maintenance, or during prescribed service operations.

The bar code scanner must be categorized as a Class 2 (II) laser.

**WARNING! LASER HAZARD.** Class 2 (II) lasers can cause damage to eyes. Avoid looking into a Class 2 (II) laser beam or pointing a Class 2 (II) laser beam into another person's eyes.

#### Workstation safety

Correct ergonomic configuration of your workstation can reduce or prevent effects such as fatigue, pain, and strain. Minimize or eliminate these effects by configuring your workstation to promote neutral or relaxed working positions.

CAUTION! MUSCULOSKELETAL AND REPETITIVE MOTION

→ HAZARD. These hazards are caused by potential risk factors that include but are not limited to repetitive motion, awkward posture, forceful exertion, holding static unhealthy positions, contact pressure, and other workstation environmental factors.

To minimize musculoskeletal and repetitive motion risks:

- Use equipment that comfortably supports you in neutral working positions and allows adequate accessibility to the keyboard, monitor, and mouse.
- Position the keyboard, mouse, and monitor to promote relaxed body and head postures.

#### Safety and electromagnetic compatibility (EMC) standards

This section provides information on:

- "U.S. and Canadian safety standards" on page 322
- "Canadian EMC standard" on page 323
- "European safety and EMC standards" on page 323
- "Australian EMC Standards" on page 328

U.S. and Canadian safety standards



The 3500 or 3500xL analyzer has been tested to and complies with standard: UL 61010-1/CSA C22.2 No. 61010-1, "Safety Requirements for Electrical Equipment for Managurament Control and Laboratory Lisa Part 1: Concrel

Equipment for Measurement, Control, and Laboratory Use, Part 1: General Requirements."

UL 61010-2-010, "Particular Requirements for Laboratory Equipment for the Heating of Materials."

The 3500 or 3500xL analyzer has been tested to and complies with the "21 CFR, 1040.10 and 1040.11 except for deviations pursuant to Laser Notice No.50, dated June 24, 2007, as applicable."

## For the Reader/Writer unit in the Applied Biosystems 3500/3500xL Genetic Analyzers

#### FCC WARNING

This device complies with Part 15 of FCC Rules. Operation is subject to the following two conditions:

- 1. This device may not cause interference, and
- 2. This device must accept any interference, including interference that may cause undesired operation of this device.

Changes or modifications not expressly approved by the party responsible for compliance could void the user's authority to operate the equipment.

#### NOTICE

This equipment has been tested and found to comply with the limits for a Class B digital device, pursuant to part 15 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference in a residential installation.

This equipment generates uses and can radiate radio frequency energy and, if not installed and used in accordance with the instructions, may cause harmful interference to radio communications. However, there is no guarantee that interference will not occur in a particular installation. If this equipment does cause harmful interference to radio or television reception, which can be determined by turning the equipment off and on, the user is encouraged to try to correct the interference by one or more of the following measures:

• Reorient or relocate the receiving antenna

- Increase the separation between the equipment and receiver.
- Connect the equipment into an outlet on a circuit different from that to which the receiver is connected.
- Consult the dealer or an experienced radio/TV technician for help.

# Canadian EMC<br/>standardThis instrument has been tested to and complies with ICES-001, Issue 3: "Industrial,<br/>Scientific, and Medical Radio Frequency Generators." Cet appareil numerique de la<br/>classe B est conforme a la norme NMB-001 du Canada.

#### Canadian Department of Communications Industry Canada (IC) Notice

This device complies with RSS-Gen of IC Rules. Operation is subject to the following two conditions:

- 1. This device may not cause interference, and
- 2. This device must accept any interference, including interference that may cause undesired operation of this device.

European safety and EMC standards EN 610

This instrument meets European requirements for safety (Low Voltage Directive 2006/95/EC). This instrument has been tested to and complies with standards EN 61010-1:2001, "Safety Requirements for Electrical Equipment for Measurement, Control and Laboratory Use, Part 1: General Requirements."

EN 61010-2-010, "Particular Requirements for Laboratory Equipment for the Heating of Materials."

EN 61010-2-081, "Particular Requirements for Automatic and Semi-Automatic Laboratory Equipment for Analysis and Other Purposes."

EN 60825-1, "Radiation Safety of Laser Products, Equipment Classification, Requirements, and User's Guide.

#### EMC

The 3500 or 3500xL analyzer meets European requirements for emission and immunity (EMC Directive 2004/108/EC).

EN 61326-1:2006 "Electrical equipment for measurement, control and laboratory use- Part 1 General EMC requirements." (Group 1, Class B)

For the Reader/Writer unit in the Applied Biosystems 3500/3500xL Genetic Analyzers

#### **CE Notice (European Union)**

Marking by the symbol indicates compliance of this ASI4000-98-BS1 RFID R/W Module to the Electromagnetic Compatibility Directive and the Low Voltage Directive of the European Union. Such making is indicative that this RFID R/W Module meets the following technical standards:

• EN 300330 – "Electromagnetic compatibility and Radio spectrum Matters (ERM); Short Range Devices (SRD)."

- EN 301489 "Electromagnetic compatibility and Radio spectrum Matters (ERM); ElectroMagnetic Compatibility (EMC) standard for radio equipment and services."
- EN 60950 "Safety of Information Technology Equipment."

Europe – CE declaration of conformity (Reader/Writer) EN 300 330-1 V1.5.1 (2006-04), EN 300 330-2 V1.3.1 (2006-04), EN 301 489-3 V1.4.1 (2002-08), EN 301 489-1 V1.6.1 (2005-09), EN 60950-1:2006

English Hereby, ART Technology Co., Ltd. declares that this ASI4000-98-BS1 is in compliance with the essential requirements and other relevant provisions of Directive 1999/5/EC.

Français

[French]

Par la présente ART Technology Co., Ltd. déclare que l'appareil

ASI4000-98-BS1 est conforme aux exigences essentielles et aux autres

dispositions pertinentes de la directive 1999/5/CE.

Deutsch

[German]

Hiermit erklärt ART Technology Co., Ltd. dass sich das Gerät ASI4000-98-BS1 in Übereinstimmung mit den grundlegenden Anforderungen und den übrigen

einschlägigen Bestimmungen der Richtlinie 1999/5/EG befindet.

Italiano

[Italian]

Con la presente ART Technology Co., Ltd. dichiara che questo

ASI4000-98-BS1 è conforme ai requisiti essenziali ed alle altre disposizioni pertinenti stabilite dalla direttiva 1999/5/CE.

Español

[Spanish]

Por medio de la presente ART Technology Co., Ltd. declara que el

ASI4000-98-BS1 cumple con los requisitos esenciales y cualesquiera otras

disposiciones aplicables o exigibles de la Directiva 1999/5/CE.

Português

[Portuguese]

ART Technology Co., Ltd. declara que este ASI4000-98-BS1 está conforme com os requisitos essenciais e outras disposições da Directiva 1999/5/CE.

Suomi

[Finnish]

ART Technology Co., Ltd. Vakuuttaa täten että ASI4000-98-BS1 tyyppinen laite on direktiivin 1999/5/EY oleellisten vaatimusten ja sitä koskevien direktiivin muiden ehtojen mukainen.

Nederlands

[Dutch]

Hierbij verklaart ART Technology Co., Ltd. dat het toestel ASI4000-98-BS1 in overeenstemming is met de essentiële eisen en de andere relevante

Česky [Czech] ART Technology Co., Ltd. tímto prohlašuje, že tento ASI4000-98-BS1 je ve shodě se základními požadavky a dalšími příslušnými ustanoveními směrnice 1999/5/ES. Dansk [Danish] Undertegnede ART Technology Co., Ltd. erklærer herved, at følgende udstyr ASI4000-98-BS1 overholder de væsentlige krav og øvrige relevante krav i direktiv 1999/5/EF. Eesti [Estonian] Käesolevaga kinnitab ART Technology Co., Ltd. seadme ASI4000-98-BS1 vastavust direktiivi 1999/5/EÜ põhinõuetele ja nimetatud direktiivist tulenevatele teistele asjakohastele sätetele. Ελληνική [Greek] ME THN ΠΑΡΟΥΣΑΑRT Technology Co., Ltd. ΔΗΛΩΝΕΙ ΟΤΙ ASI4000-98-BS1 ΣΥΜΜΟΡΦΩΝΕΤΑΙ ΠΡΟΣ ΤΙΣ ΟΥΣΙΩΔΕΙΣ ΑΠΑΙΤΗΣΕΙΣ ΚΑΙ ΤΙΣ ΛΟΙΠΕΣ ΣΧΕΤΙΚΕΣ ΔΙΑΤΑΞΕΙΣ ΤΗΣ ΟΔΗΠΑΣ 1999/5/ΕΚ. Latviski [Latvian] Ar šo ART Technology Co., Ltd. deklarē, ka ASI4000-98-BS1 atbilst Direktīvas 1999/5/EK būtiskajām prasībām un citiem ar to saistītajiem noteikumiem. Lietuvių [Lithuanian] Šiuo ART Technology Co., Ltd. deklaruoja, kad šis ASI4000-98-BS1 atitinka esminius reikalavimus ir kitas 1999/5/EB Direktyvos nuostatas. Malti [Maltese] Hawnhekk, ART Technology Co., Ltd. jiddikjara li dan ASI4000-98-BS1 jikkonforma malfid-Dirrettiva 1999/5/EC.

Magyar

[Hungarian]

Alulírott, ART Technology Co., Ltd. nyilatkozom, hogy a ASI4000-98-BS1 megfelel a vonatkozó alapvető követelményeknek és az 1999/5/EC irányelv egyéb előírásainak.

Polski

[Polish]

Niniejszym ART Technology Co., Ltd. oświadcza, że AS14000-98-BS1 jest zgodny z zasadniczymi wymogami oraz pozostałymi stosownymi postanowieniami Dyrektywy 1999/5/EC.

Slovensko

[Slovenian]

ART Technology Co., Ltd. izjavlja, da je ta ASI4000-98-BS1 v skladu z

bistvenimi zahtevami in ostalimi relevantnimi določili direktive 1999/5/ES.

Slovensky

[Slovak]

ART Technology Co., Ltd. týmto vyhlasuje, že ASI4000-98-BS1 spĺňa základné požiadavky a všetky príslušné ustanovenia Smernice 1999/5/ES.

Svenska

[Swedish]

Härmed intygar ART Technology Co., Ltd. att denna ASI4000-98-BS1 står I överensstämmelse med de väsentliga egenskapskrav och övriga relevanta bestämmelser som framgår av direktiv 1999/5/EG.

Íslenska

[Icelandic]

Hér með lýsir ART Technology Co., Ltd. yfir því að ASI4000-98-BS1 er í samræmi við grunnkröfur og aðrar kröfur, sem gerðar eru í tilskipun 1999/5/EC. Norsk

[Norwegian]

ART Technology Co., Ltd. erklærer herved at utstyret ASI4000-98-BS1 er i samsvar med de grunnleggende krav og øvrige relevante krav i direktiv 1999/5/EF.



Australian EMC Standards



This instrument has been tested to and complies with standard AS/NZS 2064, "Limits and Methods Measurement of Electromagnetic Disturbance Characteristics of Industrial, Scientific, and Medical (ISM) Radio-frequency Equipment."

## **Chemical safety**

#### General chemical safety

Chemical hazard warning

**WARNING! CHEMICAL HAZARD.** Before handling any chemicals, refer to the Material Safety Data Sheet (MSDS) provided by the manufacturer, and observe all relevant precautions.



**WARNING! CHEMICAL HAZARD.** All chemicals in the instrument, including liquid in the lines, are potentially hazardous. Always determine what chemicals have been used in the instrument before changing reagents or instrument components. Wear appropriate eyewear, protective clothing, and gloves when working on the instrument.

## Chemical safety guidelines

To minimize the hazards of chemicals:

- Read and understand the Material Safety Data Sheets (MSDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. (See "About MSDSs" on page 329.)
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing). For additional safety guidelines, consult the MSDS.
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood). For additional safety guidelines, consult the MSDS.
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended in the MSDS.
- Comply with all local, state/provincial, or national laws and regulations related to chemical storage, handling, and disposal.

#### **MSDSs**

About MSDSs Chemical manufacturers supply current Material Safety Data Sheets (MSDSs) with shipments of hazardous chemicals to new customers. They also provide MSDSs with the first shipment of a hazardous chemical to a customer after an MSDS has been updated. MSDSs provide the safety information you need to store, handle, transport, and dispose of the chemicals safely.

Each time you receive a new MSDS packaged with a hazardous chemical, be sure to replace the appropriate MSDS in your files.

Obtaining<br/>MSDSsThe MSDS for any chemical supplied by Applied Biosystems is available to you free<br/>24 hours a day. To obtain MSDSs:

- 1. Go to www.appliedbiosystems.com, click Support, then select MSDS.
- 2. In the Keyword Search field, enter the chemical name, product name, MSDS part number, or other information that appears in the MSDS of interest. Select the language of your choice, then click **Search**.
- 3. Find the document of interest, right-click the document title, then select any of the following:
  - **Open** To view the document
  - **Print Target** To print the document
  - Save Target As To download a PDF version of the document to a destination that you choose

**Note:** For the MSDSs of chemicals not distributed by Applied Biosystems, contact the chemical manufacturer.

#### Chemical waste safety

Chemical waste hazards



**CAUTION! HAZARDOUS WASTE.** Refer to Material Safety Data Sheets and local regulations for handling and disposal.

WARNING! CHEMICAL WASTE HAZARD. Wastes produced by Applied Biosystems instruments are potentially hazardous and can cause injury, illness, or death.



**WARNING! CHEMICAL STORAGE HAZARD.** Never collect or store waste in a glass container because of the risk of breaking or shattering. Reagent and waste bottles can crack and leak. Each waste bottle should be secured in a low-density polyethylene safety container with the cover fastened and the handles locked in the upright position. Wear appropriate eyewear, clothing, and gloves when handling reagent and waste bottles.

## Chemical waste safety guidelines

To minimize the hazards of chemical waste:

- Read and understand the Material Safety Data Sheets (MSDSs) provided by the manufacturers of the chemicals in the waste container before you store, handle, or dispose of chemical waste.
- Provide 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.)
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing). For additional safety guidelines, consult the MSDS.
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood). For additional safety guidelines, consult the MSDS.
- Handle chemical wastes in a fume hood.
- After emptying a waste container, seal it with the cap provided.
- Dispose of the contents of the waste tray and waste bottle in accordance with good laboratory practices and local, state/provincial, or national environmental and health regulations.
- **Waste disposal** If potentially hazardous waste is generated when you operate the instrument, you must:
  - Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
  - Ensure the health and safety of all personnel in your laboratory.

• Ensure that the instrument 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.

#### **Biological hazard safety**

## General biohazard

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. Follow all applicable local, state/provincial, and/or national regulations. Wear appropriate protective equipment, which includes but is not limited to: protective eyewear, face shield, clothing/lab coat, and gloves. All work should be conducted in properly equipped facilities using the appropriate safety equipment (for example, physical containment devices). Individuals should be trained according to applicable regulatory and company/institution requirements before working with potentially infectious materials. Read and follow the applicable guidelines and/or regulatory requirements in the following:

- U.S. Department of Health and Human Services guidelines published in *Biosafety in Microbiological and Biomedical Laboratories* (stock no. 017-040-00547-4; bmbl.od.nih.gov)
- Occupational Safety and Health Standards, Bloodborne Pathogens (29 CFR§1910.1030; www.access.gpo.gov/ nara/cfr/waisidx\_01/29cfr1910a\_01.html).
- Your company's/institution's Biosafety Program protocols for working with/handling potentially infectious materials.

Additional information about biohazard guidelines is available at: www.cdc.gov

### Safety alerts

For the definitions of the alert words **IMPORTANT**, **CAUTION**, **WARNING**, and **DANGER**, see "Safety alert words" on page xiii.

#### **Chemical alerts**

For the definitions of the alert words **IMPORTANT**, **CAUTION**, **WARNING**, and **DANGER**, see "Safety alert words" on page xiii.

General alerts for all chemicals

**WARNING!** Wear appropriate protection, including gloves, laboratory goggles, and coat whenever you work with the fluids used on this instrument, or parts that may come into contact with these fluids.

**IMPORTANT!** Use the cleaning agents as described in this manual, only. Use of cleaning agents not described in this manual can impair the instrument. Please contact your local Life Technologies sales office if you have any questions.

- Read the MSDS for this product, and follow the handling instructions.
- Avoid inhalation, contact with eyes, skin, clothing, and prolonged or repeated exposure.
- Consumables have a limited lifetime. Overusing the parts might result in poor quality data.

Specific chemical alerts

WARNING! CHEMICAL HAZARD. POP-4<sup>™</sup> POLYMER. Causes eye, skin, and respiratory tract irritation. Avoid breathing vapor. Use with adequate ventilation.

WARNING! CHEMICAL HAZARD. POP-6<sup>™</sup> POLYMER. Causes eye, skin, and respiratory tract irritation. Avoid breathing vapor. Use with adequate ventilation.



WARNING! CHEMICAL HAZARD. POP-7<sup>™</sup> POLYMER. Harmful by inhalation and if swallowed. Causes eye, skin, and respiratory tract irritation. Do NOT taste or swallow. Avoid breathing vapor (or dust). Keep container tightly closed. Use only with adequate ventilation. Wash thoroughly after handling.



**WARNING!** CHEMICAL HAZARD. Hi-Di<sup>™</sup> Formamide. Causes eye, skin, and respiratory tract irritation. Possible developmental and birth defect hazard. Avoid breathing vapor. Use with adequate ventilation.



**WARNING!** CHEMICAL HAZARD. Anode Buffer Container (ABC). May cause eye, skin and respiratory tract irritation. Avoid breathing vapor. Use with adequate ventilation.



**WARNING!** CHEMICAL HAZARD. Cathode Buffer Container (CBC). May cause eye, skin and respiratory tract irritation. Avoid breathing vapor. Use with adequate ventilation.



**WARNING!** CHEMICAL HAZARD. 1× GA Buffer/EDTA. May cause eye, skin and respiratory tract irritation. Avoid breathing vapor. Use with adequate ventilation.

#### Instrumentation alerts

General instrumentation alerts



**WARNING!** Wear appropriate protection, including gloves, laboratory goggles, and coat whenever you work with the fluids used on this instrument, or parts that may come into contact with these fluids.

Specific instrumentation alerts

fic ∠ on ∠

**WARNING!** Wear appropriate protection, including gloves, laboratory goggles, and coat whenever you work with the fluids used on this instrument, or parts that may come into contact with these fluids.

The instrument uses a Solid-state laser. Under normal operating conditions, the instrument is categorized as a Class I laser / LED product. When safety interlocks are disabled during certain servicing procedures, the laser can cause permanent eye damage, and, therefore, is classified under those conditions as a Class 3B laser.

The instrument has been tested to and complies with 21 CFR, 1040.10 and 1040.11 except for deviations pursuant to Laser Notice No.50, dated June 24, 2007, as applicable.

The instrument has been tested and complies with standard EN60825-1: 2001, "Radiation Safety of Laser Products, Equipment Classification, Requirements, and User's Guide.'



**CAUTION!** Use of controls or adjustments or performance of procedures other than those specified herein may result in hazardous radiation exposure.

#### Laser Parameter

Wave length 505nm, Output power 20mW

#### **LED** Parameter

Emitting color Natural White, Luminous Intensity 250 Cd.

## Documentation

## **Related documentation**

The following related documents are shipped with the system:

Document	Part number	Description	
Applied Biosystems 3500/3500xL Genetic Analyzers Quick Reference Card	4401662	<ul> <li>Provides a flowchart on how to run your samples and instrument</li> <li>Provides a table of maintenance tasks and</li> <li>Contains Data Collection Software reference guide</li> </ul>	
Applied Biosystems 3500 Series Genetic Analyzer Site Preparation Guide (4401689)	4401663	Provides information about the space, environmental, and electrical requirements needed to support the Applied Biosystems 3500/3500xL Genetic Analyzers.	
<b>Note:</b> The purpose of the Site Prep Guide is to help you prepare your site for installation of the 3500 or 3500xL analyzer. For specific details about your system, please refer to this user guide.			

Portable document format (PDF) versions of this guide as well as the Quick Reference Card and the Warranty statement are also available on the Applied Biosystems 3500/3500xL Genetic Analyzers the software installation CD, which will be shipped with the system.

Note: For additional documentation, see "How to obtain support" on page xvii.

### Obtaining information from the Help system

The 3500 Series Data Collection Software interface has instructions guiding the user through basic tasks of the workflow and expanded help information for complex decisions and operations. Users can access these instructions by clicking the help icon ?

The 3500 or 3500xL analyzer has a Help system that describes how to use each feature of the user interface. Access the Help system by doing one of the following:

- Click **?** in the screens of the 3500 Series Data Collection Software window.
- Select Help > Help Contents.

You can use the Help system to find topics of interest by:

- Reviewing the contents
- Searching for a specific topic
- Searching an alphabetized index

### Send us your comments

Applied Biosystems welcomes your comments and suggestions for improving its user documents. You can e-mail your comments to:

techpubs@appliedbiosystems.com

**IMPORTANT!** The e-mail address above is for submitting comments and suggestions relating *only* to documentation. To order documents, download PDF files, or for help with a technical question, see "How to obtain support" on page xvii.

## Index

#### Symbols

```
.annotation.txt 36, 87
.csv 75
.fsta 36, 87
.html 87
.pdf 87
   action log 209
   audit reports 214
   e-sig report 222
   fragment/HID install performance check 138
   plate layout 50
   sample quality reports 95
   sequencing install performance check 128
   spatial calibration report 102
   spectral calibration 116
   trace quality reports 85
   user report 206
.phd.1 36, 87
.qual 36, 87
.scf 36, 87
.seq 36, 87
.txt 75, 87
.xls 75, 87
.xml 75
```

#### **Numerics**

3500 Series Data Collection Software 3500 Software 25 compatible secondary analysis applications 14 Daemon 24 Dashboard 15 default settings 32, 33 files generated 14 Library workflow 17  $\log in 26$ Main workflow 16 Maintenance workflow 17 navigate 19 overview 14 parts of 15 required procedure to start 24 Server Monitor 25 status icon in Windows tray 24

use without an instrument 19 8-tube strip plate base 53 plate type 44, 51, 105, 120, 145 96 Fast tube, plate type 145

#### Α

ABC. See anode buffer abort injection 69 account setup 199 suspended, activate 203 suspension 199, 226 user 201 Acronyms xvi action log contents 212 display 209 Administrator user role 203 allelic ladder internal validation 52 re-injection 67 requirements for accurate genotyping 52, 68 results group for one allelic ladder per run folder 52 run requirements 155 sample name for re-injection 49 sample type 48 amber indicator light 23 amplicon, specifying 49 Analysis Range fragment analysis 181 HID analysis 186 anode buffer change 237 described 9 on-instrument limits 30 part number 9 storage conditions 9 troubleshoot 303 AnyDye calibrate spectral 108 create dye set 169 Applied Biosystems customer feedback on documentation 336

Information Development department 336 archive audit records 214 data files 255 library items 254 array See also capillary array view 77 array selection, assign plate contents 71 as-needed instrument maintenance 232 assavs assign to plate 49, 73 create 147 defined 46, 147 instrument protocol 150 normalize data 49 assign plate contents amplicon and specimen 49 array selection (capillary-to-plate map) 71 assay 49, 73 file name convention 49, 73 name samples 48, 70 overview 46 parts of screen 46 plate view 48, 70 plate view, customize 71 results group 49, 73 sample type 48table view 71 template for 75 Assume Linearity 183 audit records factory-provided items, creation date 142 view in library 142 audit, administrators action log 209, 212 archive records 214 audit actions 210, 212 audit mode 208 audit reason settings 208 audit records, object 210 audited objects and actions 208 enable or disable 207 export records 215 export settings 224 import settings 224 object audit history 209, 210 overview 197 purge records 214 restore records 214 system configuration history 209, 211 troubleshoot 312 when security is disabled 207

audit, users enter reason for change 227 overview 225 view audit records 142 auto-analysis plate setting 146 average quality value (QV), monitor run 64

#### В

barcode 145 barcode scanner 8 basecall probability of error 84 Basecaller 175 basecaller version in basecalling protocol and View Sequence results 80 basecalling protocols create 174 defined 174 export 141 import 141 in assay 150 Basepair Accuracy 125 BigDye® Terminator kits and standards 259 biohazardous waste, handling 331 bold text, when to use xvi borrowing defined 112 limits 34 Broad Peak (BD) and SQ 188 flag in view fragment/HID results 89 threshold 188 buffer, anode. See anode buffer buffer, cathode. See cathode buffer Buffer-Pin Valve 28

#### С

calibration See spatial calibration See spectral calibration capillary array change 252 check stored 240 described 12 fill with polymer 251 on-instrument limits 30 part number 13 storage conditions 240 capillary-to-plate map diagram 51 display in software 71 cathode buffer change 238 described 9 on-instrument limits 30 part number 9 storage conditions 9 troubleshoot 304 CAUTION, description xiii CBC. See cathode buffer Change Polymer Type maintenance wizard 247 check valve 28 chemical safety 328 chemical waste safety 330 Clear range 176 computer maintenance. See maintenance, computer name requirement 22 start up 24 condition number, spectral calibration 109 conditioning reagent described 11 part number 11 place on instrument 250 consumables anode buffer 9, 237 capillary array 240, 252 cathode buffer 9, 238 conditioning reagent 11, 250 fragment analysis and HID 260 Hi-Di<sup>™</sup> Formamide 12 on-instrument limits 30 polymer 10, 245, 247, 249 sequencing analysis 259 status in Dashboard 29, 236 contiguous read length. See CRL controls, specifying 48 conventions bold text xvi for describing menu commands xvi **IMPORTANTS!** xvi italic text xvi Notes xvi user attention words xvi CRL ambient temperature dependence 30 defined 81 distribution report 85 performance check 125 QV settings 178 report 85 result 81 threshold 178

CRL Basepair Accuracy 125 CRL Median and SD 125 custom dye set. *See* dye sets, custom customer feedback, on Applied Biosystems documents 336 customize sample info 48

#### D

daily instrument maintenance 230, 231 DANGER, description xiii Dashboard 29 access from Main workflow 19 consumables status 236 instrument status 53 troubleshoot 309 data collection software. See 3500 Series Data Collection Software data files. See sample data files data troubleshooting 306 datastore back up during uninstall 254 date format, setting 33 default settings, specifying 32, 33 define plate properties 44 detection cell temperature 42, 104, 119, 129 documentation, related 335 duplicate injection See also re-injection defined 60preview run 60duplicate library item 140 dye set, overview.See dye sets dye sets calibrate 103 create 168 custom, calibrate 108 custom, create 169 defined 168 export 141 fragment analysis 261 HID analysis 262 import 141 sequencing 261

#### Ε

electrical safety 319 electromagnetic compatibility standards. *See* EMC standards electronic signature records actions 221 view in library 142 electronic signature, administrators

actions that allow e-sig 217 enable or disable 216 export 223 export settings 224 functions that require e-sig 217 import settings 224 is signed field 228 overview 197 report 222 when security is disabled 216 electronic signature, users is signed field 228 overview 225 signing 227 electropherogram troubleshooting 306 EMC standards 322 EPT view 78 ergonomics, safety 321 error messages display detail 312 instrument 299, 313 link plate 304 RFID 304 spatial calibration 300 spectral calibration 301 e-sig. See electronic signature export assays 141 audit records 215 audit settings 224 basecalling protocols 141 dye sets 141 e-sig records 223 e-sig settings 224 file name convention 141 fragment analysis protocols 141 HID analysis protocols 141 instrument protocols 141 plate 75, 141 protocol 141 QC protocols 141 results groups 141 security settings 224 sequencing analysis protocols 141 sequencing results, reports, traces 87 size standards 141 sizecalling protocols 141 spatial calibration 101 spectral calibration 115 user account settings 224

#### F

factory-provided library items 139 fast plate plate base 53 plate type 44, 51, 105, 120, 145 file location default 35 in file name convention 154 with file name convention 50, 151with results group 50, 155 without file name convention 50, 151 without results group 50, 155file name conventions assign to plate 49, 73 create 151 defined 151 export 141 file location 154 import 141 in plate 151 file name format default format 50 with file name convention 151 without file name convention 50, 151 file name maximum length 154 fill and fill series, sample name 48, 71, 72 Fill Array with Polymer maintenance wizard 251 first run migration effects 53 flags display flag table in monitor run 62 monitor run 62 re-injecting samples with 65 Flat Profile 176 formamide, sample type for 48 fragment analysis dye sets 261 reagents, part numbers, storage conditions 260 run modules 264 fragment analysis protocols create 193 defined 193 export 141 import 141 in assay 150 fragment analysis results. See review results, fragment/HID fragment/HID install performance check analyze data in secondary analysis software 136 estimated run time 129 evaluate data 135 examples 137 historical reports 138

history 137 instrument, prepare 129 plate, prepare 129 recommended monthly schedule 231 report 137 run 132 signing 228 standards 130 troubleshoot 136, 303 what occurs 134 when to perform 129

#### G

GeneMapper ID-X protocols See also HID analysis protocols create 195 GeneMapper protocols See also fragment analysis protocols create 193 green indicator light 23 GS600 LIZ® v2.0 size standard 171, 260 GS600(60-600)LIZ+Normalization 171 GS600(80-400)LIZ+Normalization 171 GS600LIZ+Normalization 171 guidelines chemical safety 328 chemical waste disposal 330 chemical waste safety 330

#### Η

hazard icons. See safety symbols, on instruments 315 hazard symbols. See safety symbols, on instruments hazards. See safety Help system, accessing 336 HID analysis allelic ladder validation 52 dye sets 262 reagents, part numbers, storage conditions 260 results group for one allelic ladder per run folder 52 results. See review results, fragment/HID run modules 264 HID analysis protocols create 195 defined 195 export 141 import 141 in assay 150 HID install performance check. See fragment/HID install performance check Hi-Di<sup>™</sup> Formamide on-instrument limits 12

part number 12 Hi-Di<sup>™</sup> sample type 48

#### I

import assays 141 audit settings 224 basecalling protocols 141 dye sets 141 e-sig settings 224 file name convention 141 fragment analysis protocols 141 HID analysis protocols 141 instrument protocols 141 plate 44, 75, 141 QC protocols 141 results group 141 sample files 81 sample files, fragment/HID 89 security settings 224 sequencing analysis protocols 141 size standards 141 sizecalling protocols 141 user account settings 224 IMPORTANT, description xiii Information Development department, contacting 336 injection defined 57 duplicate 60 folder 159 re-injection 65 replicate 60 specify new instrument protocol 65 injection list abort injection 69 blank 60 columns in table 62completed injections 62 duplicate injection 60 modify 60 monitor run 61 preview run 59 re-injection 65 replicate injection 60 samples with assays that specify more than one instrument protocol 61 terminate 69 Install Capillary Array maintenance wizard 252 installation category 319 instrument error messages 313 load plate 53 maintenance. See maintenance

move and level 243 name, specifying 33 prepare for run 42 reset 314 routine cleaning 242 shutdown 253 start up 22 status indicator lights 23 status, in Dashboard 53 theory of operation 5troubleshoot 299, 313 verification. See fragment/HID performance check, sequencing performance check instrument operation 236 instrument operation, safety 318 instrument protocols create 165 defined 165 export 141 import 141 in assay 150 normalization settings 167 instrument run abort injection 69 defined 57 monitor 61 pause 69 preview 61 resume 69 start 61, 69 start with existing plate 55 stop 69 terminate injection list 69 instrument run name See also run name folder name 159 instrument run views 77 instrument sensor details view 305Instrument Shutdown maintenance wizard 253 IQ/OQ runs. See fragment/HID performance check, sequencing performance check is signed field 228 italic text, when to use xvi

#### L

label peaks 93, 94 laser classification 320, 321 laser safety 320 bar code scanner 321 requirements 320, 321 libraries

assays 147 basecalling protocols 174 dye sets 168 file name conventions 151 fragment analysis protocols 193 HID analysis protocols 195 instrument protocols 165 MicroSeq ID analysis protocols 191 QC protocols 184 results group 156 sequencing analysis protocols 189 size standards 171 sizecalling protocols 179 library AB 139 access 140 archive 254 create new item 140 delete entry 141 delete entry, auditing 141 edit entry 141 export entry 141 factory-provided 139 factory-provided, create new item from 140 file name conventions 151 filter 142 GeneMapper ID-X protocols 195 import entry 141 locked 139 locked, create new item from 140 overview 196 search 142 SeqScape protocols 189 sort 142 template 139 template, create new item from 140 view audit records 142 view e-sig records 142 linearity, large-fragment size standards 183 link plate 54, 57 link plate error message 304 link plate troubleshooting 309 load plate for run 56 load plate troubleshooting 309 locked library item 139 log files, search and use 305 log in 26, 225 logged-in user name determining 226 display 203 in user account 202

#### Μ

Main workflow access from Dashboard 19 maintenance calendar create entries 233 default entries 233 F entries 233 FR entries 233 recommended entries 233 view 232 maintenance notifications complete a task 28, 229 Dashboard 28, 229 dismiss a task 28, 229 log 234, 257 trigger time, setting 33 maintenance wizards change polymer type 247 fill array with polymer 251 install capillary array 252 instrument shutdown 253 overview 244 remove bubbles 251 replenish polymer 245 wash pump and channels 249 maintenance, computer data files, archive 255 defragment 257 disk space, monitor 256 library items, archive, purge, and restore 254 uninstall software 254 maintenance, consumables anode buffer, change 237 capillary array, change 252 capillary array, check stored 240 cathode buffer, change 238 polymer, change type 247 polymer, fill array 251 polymer, replenish 245 polymer, store partially used pouch 249 maintenance, instrument annual 232 as-needed 232 daily 230 monthly 231 move and level the instrument 243 planned 232 pump, flush water trap 241 pump, remove bubbles 251 pump, wash chamber and channels 249 quarterly 231 routine cleaning 242 schedule 229

shutdown 253 weekly 231 workflow 17 manual commands troubleshoot 313 matrix standard part numbers 260 menu commands, conventions for describing xvi metric analysis display results 81 flag settings 177 warnings 81 MicroSeq ID analysis protocols create 191 defined 191 migration effects, first run 53 Min. Peak Half Width 182, 187 mixed base color 37 QV range 37 settings 84 threshold 176 Mobility file 175 monitor run current run 61 flag settings 177 flags in injection list 62 flags in plate view 62injection list 61 injection list, modify 61 re-injections 65 start, stop, resume run 69 troubleshoot 311 monthly instrument maintenance 231 move the instrument 243 moving and lifting safety computers and monitors 317 instrument 317 moving parts, safety 318 MSDSs about xiii description 329 obtaining xvii, 329

#### Ν

negative control 48 non-linearity in large fragment size standards 183 normalization and SQ (sizing quality) 183, 188 factor 91, 167 factor in secondary analysis software 91 factor thresholds 91, 167 flag in view fragment/HID results 89 GS600 LIZ® v2.0 size standard 171 how the software normalizes 91 overview 7 review data 90 size standards 171 target 91, 167 notifications log for maintenance actions 234, 257 maintenance in Dashboard 28, 229 SAE module 200

#### 0

object audit history contents 210 display 209 offscale flag monitor run 64 view fragment/HID results 89 online Help. *See* Help system oven temperature in instrument protocol 167 pre-heat in Dashboard 42 overlay report 95 overvoltage category (rating) 319 owner, plate 145

#### Ρ

part numbers anode buffer 9 capillary array 13 cathode buffer 9 conditioning reagent 11 fragment/HID analysis standards 260 Hi-Di<sup>™</sup> Formamide 12 sequencing analysis reagents 259 password change your own 226 expiration 199 restrictions 199 pause instrument run 69 Peak Amplitude Thresholds 181, 187 Peak Window Size 182 pending plates 76 performance check See fragment/HID install performance check See sequencing install performance check permissions, user account 203, 225 physical hazard safety 318 planned maintenance 232 plate See also assign plate contents

assay 49, 73 assembly 53 base, for standard versus fast and 8-strip 53 create 43 create for import 73 create from template 43 edit 74 edit existing 46 export 75 file name convention 49,73import 44, 75 link 54, 57 load in instrument 53 load in software 56 overview 143 owner 145 pending 76 print 50 processed 76 properties 44 results group 49, 73 sample plate, prepare 51 select for run 54 setup error 309signing 228 type 145 unlink 57 plate base, standard versus fast and 8-tube strip 44, 51, 53, 105, 120, 145 plate definition 143 plate grid report 50 plate import template 73 plate layout .pdf 50 print 50 saving for future use 75 plate map, print 50 plate properties 145 plate record. See plate plate report fragment/HID 95 sequencing, generate 85 plate report, sequencing quality ranges 38 plate template create 75 create plate from 43 save well attributes 75 plate type default 34 set default 76 specify 44, 51, 105, 120, 145 plate view

assign plate contents 48 customize assign plate contents 71 monitor run 62 plate-to-capillary map diagram 51 display in software 71 plots, fragment/HID display plots 91 label peaks 93 overlay 93 scale 92 settings 92 settings, default 35 zoom 92 polymer ambient temperature requirements 30 change type 247 described 10 expiration 11 fill array 251 on-instrument limits 30 oven temperature, POP-4, POP-6, POP-7 42 part numbers 11 replenish 245 storage conditions 10 storing partially used pouch 249 Polynomial Degree 182 POP polymer. See polymer positive control 48 preferences borrowing event limits 34 date format 33 file location 35instrument name 33 maintenance notification time 33 overview 32plate type 34 plot settings 35 reports 34 reports, sequencing 38 sequencing export 36sequencing trace 36system 32 table settings 35 user 33 pre-heat duration 42 POP-4, POP-6, POP-7 polymer 42 pressurized fluids, safety 318 preview run duplicate injection 60 injection list 59 start instrument run 61

primer peak, eliminating 181 print action logs 209 audit reports 213 e-sig report 222 fragment/HID install performance check 137 plate layout 50 sample quality reports 95 sequencing install performance check 128 spatial calibration report 102 spectral calibration 116 trace quality reports 85 user report 205 processed plates 76 Pull-Up peak flag 89 settings 183 pump flush water trap 241 remove bubbles 251 wash chamber and channels 249 PUP Score 81 pure base color 37 QV range 37 pure base settings 84 purge audit records 214 library items 254

#### Q

QC protocols create 184 defined 184 export 141 import 141 in assay 150 QC report, sequencing generate 85 quality ranges 38 quality flags fragment/HID 62, 89 sequencing 62, 81 quality settings fragment analysis 183 HID analysis 188 sequencing analysis 177 quality value See also QV probability of error 84 spectral calibration 109 quarterly instrument maintenance 231 Quick Start 55 QV (average quality value) See also average quality value monitor run 64 QV settings, sequence basecalling protocol 177 color and ranges 37 monitor run screen 177 recommended ranges 84 QV20+ report 85 result 81 setting 178

#### R

radioactive waste, handling 331 Read Length 125 recent plates 76 runs 76 red indicator light 23 re-injection See also duplicate injection allelic ladder 67 collect data for specific wells 65 defined 65 file location 159 in view sequencing results 85 instrument protocol, selecting 65 monitor run 65 order 65 samples, all in injection 65 samples, individual 65 re-injection button graved 65 re-injections in view fragment/HID results 95 Remove Bubbles maintenance wizard 251 rename samples in review results 97 repetitive motion, safety 321 Replenish Polymer maintenance wizard 245 replicate injection. See injection list report settings fragment/HID 95 sequence 86 reports action log 209, 212 audit 213 CRL and CRL distribution 85 default settings 34 default settings, sequencing 38 e-sig 222

export sequencing 87 fragment/HID install performance check 137 fragment/HID results 95 object audit history 209, 210 plate grid 50 sample quality 95 sequencing defaults 38 sequencing install performance check 127 sequencing results 85 spatial calibration 102 spectral calibration 116 system configuration history 209, 211 trace score 85 user 205 reset instrument 314 restore audit records 214 library items 254 results See review results See review results, fragment/HID analysis See review results, sequencing results groups assign to plate 49, 73 create 156 defined 155 export 141 folder 159 for HID applications 52, 155, 161 import 141 store samples by plate name 160 resume instrument run 69 review results currently running plate, fragment/HID 88 currently running plate, sequencing 80 import sample files 81 import sample files, fragment/HID 89 previously run plate 81 previously run plate, fragment/HID 89 rename 97 show/hide columns 81 sort 81 review results, fragment/HID access 88 label peaks 93 modify data in secondary analysis software 98 normalized data 90 overlay plots 93 plots 91 quality flags 89 re-injections 95 sample quality 89 scale 92

sizing 94 sizing curve, overlay 95 thumbnails 94 troubleshoot 311 zoom 92review results, sequencing 87 See also sequence quality access 80 base colors 37export defaults 36 metric analysis 81 modify data in secondary analysis software 98 QV colors and ranges 37 re-inject 85 thumbnails 84 trace defaults 36trace quality reports 85 RFID error message 304 troubleshoot 304 run module 166 run modules fragment/HID analysis 264 sequencing analysis 263 run name default 57 enter 57 run voltage 167 run, see instrument run

## S

SAE module See audit See electronic signature See security safety bar code scanner 321 before operating the instrument 317biological hazards 331 chemical 328 chemical waste 330 electrical 319 ergonomic 321 guidelines 328, 330 instrument operation 318 laser 320, 321 moving and lifting computers and monitors 317 moving and lifting instrument 317 moving parts 318 physical hazard 318 pressurized fluids. See also compressed gases, safety repetitive motion 321

solvents 318 standards 322 workstation 321 safety labels, on instruments xiv, 317 safety standards 322 safety symbols, on instruments 315 sample data file storage file name convention location 50 injection folder 159 instrument run name folder 159 results group location 50 results group name folder 159 sample data file storage location 18 sample data files, archive 255 sample info, enter 48 sample name allelic ladder 49 assign 48, 71 fill and fill series 48, 71 sample plate. See plate sample quality See also review results, fragment/HID in monitor run 62in review results 89 sample type plate view 48 table view 71 sample view 77 scale plots 92 scientist user role 203 secondary analysis fragment 193 GeneMapper 193 GeneMapper ID-X 195 HID 195 MicroSeq ID 191 SeqScape 189 sequencing 189, 191 security policies 199 security, administrators account setup 199 disable, effect on audit and e-sig 198 enable or disable 198 export settings 224 export user account settings 224 import settings 224 import user account settings 224 notification 199, 200 overview 197 security policies 199 spaces in user names 200 user accounts 201

user report 205 user role 203username restrictions 199 security, users account suspension 226 log in 225 overview 225 password change 226 permissions 225 session timeout 226 SeqScape protocols See also sequence analysis protocols create 189 sequence quality See also review results, sequencing in monitor run 62 in review results 81 mixed base settings 84 pure base settings 84 recommended ranges 84 trace quality reports 85 sequencing analysis dye sets 261 run modules 263 sequencing analysis protocols create 189 defined 189 export 141 import 141 in assay 150 sequencing analysis reagents, part numbers and storage conditions 259 sequencing install performance check evaluate data 126 examples 127 historical reports 128 history 127 instrument, prepare 119 plate, prepare 120 recommended monthly schedule 231 report 127 required well positions 120 run 122 signing 228 standards 120 thresholds 125 troubleshoot 302 what occurs 124 when to perform 119 sequencing results. See review results, sequencing Server Monitor 25 service log 258 session timeout 199, 226

shutdown 253 signal strength report, sequencing 85 signing, electronic signature 227 Size Calling Method fragment analysis 181 HID analysis 186 size standard plot 94 size standards create 171 defined 171 export 141 import 141 in QC protocol 186 in sizecalling protocol 180 large fragment, non-linearity 183 normalization 171 part numbers 260 sizecalling protocols create 179 defined 179 export 141 import 141 in assay 150 sizing curve, overlay 95 sizing quality. See SQ (sizing quality) Sizing Range fragment analysis 181 HID analysis 186 sizing report 95 sizing results fragment/HID, label peaks 94 Slope Thresholds Peak Start and End 182, 187 Smoothing 181, 186 software without instrument 19 software. See 3500 Series Data Collection Software solvents, safety 318 sort library 142 multi-column 142 spatial calibration export 101 historical reports 102 perform 99 purpose 99 signing 228 troubleshoot 300 when to perform 99 specimen, specifying 49 spectral calibration AnyDye 108 borrowing event limits 34 capillary sharing 112 condition number 109

estimated run time 103 evaluate data 110 examples 114 export 115 historical reports 116 history 117 instrument, prepare 104 not retained for new polymer type 106 perform 106 plate, prepare 104 pull-down peaks 112 purpose 103 quality value 109 report 116 required well positions 105 signing 228 standards 104 troubleshoot 301 what occurs 112 when to perform 103zoom 111 spectral pull-up flag 89 SQ (sizing quality) and broad peaks 188 and normalization 183, 188 flag in monitor run 64 flag in view fragment/HID results 90 fragment analysis 183 fragment analysis, difference from GeneMapper ID-X 183 HID analysis 188 monitor run 64 weighting 188 stand-alone installation of software 19 standard plate plate base 53 plate type 44, 51, 105, 120, 145 standards EMC 322 fragment/HID install performance check 130 safety 322 sequencing install performance check 120 spectral calibration 104 start instrument run 69 start the system 21 status consumables in Dashboard 26 front panel indicator lights 23 software, icon in Windows tray 24 stop instrument run 69 symbols, safety 315 system configuration history contents 211

display 209 system defaults date format 18 system description 1 system start up 21

#### Т

Table of Acronyms xvi table view, assign plate contents show/hide columns 72 sort 72 use 71 user-defined fields 48 tables results, show/hide columns 81 settings 72 settings, default 35 show/hide columns 72 sort 97 sort, multi-column 72 technician user role 203 template create plate from 43 library 139 plate 75 plate import 73 terminate instrument run 69 theory of operation 5thumbnails fragment/HID 94 sequencing 84 timeout, session 199, 226 trace quality reports 85 Trace Score **81**, 178 Trace Score report 85 training, information on xvii troubleshoot anode buffer container 303 audit 312 cathode buffer container 304 Dashboard 309 data 306 electropherogram 306 fragment/HID install performance check 303 instrument 299, 313 link plate 304, 309 load plate 309 log files, search and use 305manual commands 313 monitor run 311 red indicator status light 23 review results, fragment/HID 311

RFID 304 sequencing install performance check 302 spatial calibration 300 spectral calibration 301 view instrument sensor details 305 True Profile 176

#### U

uninstall the software 254 Uninterruptible Power Supply 8 unlink plate 57 user account activate suspended 203 create or edit 201 delete 203 inactivate 203 permissions 203 user attention words, described xvi user role, create 203 user-defined fields displayed in table view only 48 including in file name 154 specifying for samples 48

#### V

validation, allelic ladder 52 verification runs. *See* fragment/HID performance check, sequencing performance check view fragment/HID results. *See* review results, fragment analysis results, HID analysis results view run results. *See* review results view sequencing results. *See* review results, sequencing results

#### W

WARNING, description xiii warnings, metric analysis 81 Wash Pump and Channels maintenance wizard 249 waste disposal, guidelines 330 waste profiles, description 330 weekly instrument maintenance 231 weighting, SQ 188 well attributes assign plate contents 46 save in template 75 workflow diagram 21, 79 workstation safety 321

#### Ζ

zoom

fragment/HID plots 92 plate view 71 spectral calibration 111

Part Number 4401661 Rev. C 06/2010



Headquarters

850 Lincoln Centre Drive | Foster City, CA 94404 USA Phone 650.638.5800 | Toll Free 800.345.5224 www.appliedbiosystems.com

#### International Sales

For our office locations please call the division headquarters or refer to our Web site at www.appliedbiosystems.com/about/offices.cfm