

thermo scientific



Thermo

Compound Discoverer User Guide

Software Version 3.1

XCALI-98120 Revision A • July 2019



© 2019 Thermo Fisher Scientific Inc. All rights reserved.

Compound Discoverer, Exactive, Foundation, FreeStyle, Mass Frontier, Metabolika, mzCloud, mzVault, Normalized Collision Energy, Orbitrap ID-X, Q Exactive, ToxID, and TraceFinder are trademarks; Orbitrap, Orbitrap Fusion, Orbitrap Lumos, Thermo Scientific, and Xcalibur are registered trademarks of Thermo Fisher Scientific Inc. in the United States.

The following are registered trademarks in the United States and other countries: Adobe, Flash, and Reader are registered trademarks of Adobe Systems Incorporated. Microsoft, Internet Explorer, Visual C++, Windows, and Excel are registered trademarks of Microsoft Corporation. Xeon is a registered trademark of Intel Corporation. ActiveReports is a registered trademark of GrapeCity, Inc.

ChemSpider is a registered trademark of ChemZoo Inc. KEGG is a registered trademark of Kanehisa, Minoru (an individual).

All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Thermo Fisher Scientific Inc. provides this document to its customers with a product purchase to use in the product operation. This document is copyright protected and any reproduction of the whole or any part of this document is strictly prohibited, except with the written authorization of Thermo Fisher Scientific Inc.

The contents of this document are subject to change without notice. All technical information in this document is for reference purposes only. System configurations and specifications in this document supersede all previous information received by the purchaser.

This document is not part of any sales contract between Thermo Fisher Scientific Inc. and a purchaser. This document shall in no way govern or modify any Terms and Conditions of Sale, which Terms and Conditions of Sale shall govern all conflicting information between the two documents.

Release history: Revision A July 2019

Software version: Compound Discoverer 3.1.x

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

Contents

	Preface
	Related Documentation xvii
	System Requirements
	Installation Instructions xxii
	Installing the Compound Discoverer Application and the mzVault
	Libraries
	Installing the mzVault 2.2 Applicationxxiv
	Special Notices
	Contacting Usxxvi
Chapter 1	Introduction
	New Features and Enhancements
	New Features
	Enhancements
	Supported File Formats
	Starting the Application
	Choosing the Toolbar Icon Size6
	The Application Window7
	Application Menu Bar7
	Application Toolbar
	Understanding Processing Workflows
	Targeted Workflows for Expected Compounds
	Untargeted Workflows for Identifying Unknown Compounds
	Marking Background Compounds in an Untargeted Analysis
	Nomenclature for the Defined Processing Workflows
	The Defined Processing Workflows
	Best Scans for Composition Prediction and Spectral Matching
	Best MS1 Scan for Isotope Pattern Matching
	Best MS2 Scan for Fragments Matching and Spectral Comparison 33
	Quality Control Samples for Batch Normalization
	Stable Isotope Labeling
	FISh Scoring for Targeted Compounds and Proposed Structures
	Calculating the Mass Defect of an Elemental Composition

C

	Using mzLogic to Score Candidates for Unknown Compounds	39
	Running an Untargeted Analysis that Includes the mzLogic Node	
	Running an mzLogic Analysis from the mzLogic Analysis View	40
	Managing the Start Page's Recent File Lists	40
	Keeping a File Name at the Top of a Recent File List	40
	Removing Files from the Recent File Lists	40
	Exploring File Folders from the Recent File Lists.	
	Auto-Hiding the Start Page, the Chromatograms View, and the Mass	
	Spectrum View	41
	Working with the Tabbed Documents	41
	Tabbed Documents List	
	Opening a Hidden Tabbed Document	
	Tab Groups	
	Shortcut Menu Commands for Tabbed Documents	
	Rearranging the Tabbed Documents and Graphical Views	
	Moving a Graphical View to Another Monitor	
	Using the Mouse Pointer to Rearrange Tabbed Documents and	
	Graphical Views	45
	Using the Shortcut Menu Commands to Rearrange the Tabbed	
	Documents.	47
Chapter 2	Getting Started	
	Setting Up the Lists and Libraries and Checking Access to the Online	
	obtime op me ziele und zielande und entering riebete to the omine	
	Databases	49
	Databases	
	Databases	50
	Databases	
	Databases Setting Up the Expected Compounds Library for a Targeted Analyses Checking the Access to the Online Databases for Identifying Unknown Compounds Setting Up a Study and an Analysis and Processing a Data Set Setting Up a New Study and a New Analysis by Using the Wizard Troubleshooting an Analysis	
	Databases Setting Up the Expected Compounds Library for a Targeted Analyses Checking the Access to the Online Databases for Identifying Unknown Compounds Setting Up a Study and an Analysis and Processing a Data Set Setting Up a New Study and a New Analysis by Using the Wizard Troubleshooting an Analysis Submitting a Run to the Job Queue	
	Databases Setting Up the Expected Compounds Library for a Targeted Analyses Checking the Access to the Online Databases for Identifying Unknown Compounds Setting Up a Study and an Analysis and Processing a Data Set Setting Up a New Study and a New Analysis by Using the Wizard Troubleshooting an Analysis Submitting a Run to the Job Queue Opening the Result File, Reviewing the Data, and Creating a Report	
	Databases Setting Up the Expected Compounds Library for a Targeted Analyses Checking the Access to the Online Databases for Identifying Unknown Compounds Setting Up a Study and an Analysis and Processing a Data Set Setting Up a New Study and a New Analysis by Using the Wizard Troubleshooting an Analysis Submitting a Run to the Job Queue Opening the Result File, Reviewing the Data, and Creating a Report Opening the Result File from the Job Queue	
	Databases Setting Up the Expected Compounds Library for a Targeted Analyses Checking the Access to the Online Databases for Identifying Unknown Compounds Setting Up a Study and an Analysis and Processing a Data Set Setting Up a New Study and a New Analysis by Using the Wizard Troubleshooting an Analysis Submitting a Run to the Job Queue Opening the Result File, Reviewing the Data, and Creating a Report Opening the Compounds that the Analysis Detected or Found.	
	Databases Setting Up the Expected Compounds Library for a Targeted Analyses Checking the Access to the Online Databases for Identifying Unknown Compounds Setting Up a Study and an Analysis and Processing a Data Set Setting Up a New Study and a New Analysis by Using the Wizard Troubleshooting an Analysis Submitting a Run to the Job Queue Opening the Result File, Reviewing the Data, and Creating a Report Opening the Result File from the Job Queue Reviewing the Compounds that the Analysis Detected or Found. Displaying Only the Items of Interest in a Result Table.	
	Databases Setting Up the Expected Compounds Library for a Targeted Analyses Checking the Access to the Online Databases for Identifying Unknown Compounds Setting Up a Study and an Analysis and Processing a Data Set Setting Up a New Study and a New Analysis by Using the Wizard Troubleshooting an Analysis Submitting a Run to the Job Queue Opening the Result File, Reviewing the Data, and Creating a Report Opening the Compounds that the Analysis Detected or Found. Displaying Only the Items of Interest in a Result Table. Printing a Report Using a Defined Template.	
Chapter 3	Databases Setting Up the Expected Compounds Library for a Targeted Analyses Checking the Access to the Online Databases for Identifying Unknown Compounds Setting Up a Study and an Analysis and Processing a Data Set Setting Up a New Study and a New Analysis by Using the Wizard Troubleshooting an Analysis Submitting a Run to the Job Queue Opening the Result File, Reviewing the Data, and Creating a Report Opening the Result File from the Job Queue Reviewing the Compounds that the Analysis Detected or Found. Displaying Only the Items of Interest in a Result Table. Printing a Report Using a Defined Template.	
Chapter 3	Databases Setting Up the Expected Compounds Library for a Targeted Analyses Checking the Access to the Online Databases for Identifying Unknown Compounds Setting Up a Study and an Analysis and Processing a Data Set Setting Up a New Study and a New Analysis by Using the Wizard Troubleshooting an Analysis Submitting a Run to the Job Queue Opening the Result File, Reviewing the Data, and Creating a Report Opening the Compounds that the Analysis Detected or Found. Displaying Only the Items of Interest in a Result Table. Exporting a List of Compounds to a Spreadsheet Printing a Report Using a Defined Template. Starting a New Study and Setting Up an Analysis	
Chapter 3	Databases Setting Up the Expected Compounds Library for a Targeted Analyses Checking the Access to the Online Databases for Identifying Unknown Compounds Setting Up a Study and an Analysis and Processing a Data Set Setting Up a New Study and a New Analysis by Using the Wizard Troubleshooting an Analysis Submitting a Run to the Job Queue Opening the Result File, Reviewing the Data, and Creating a Report Opening the Result File from the Job Queue Reviewing the Compounds that the Analysis Detected or Found. Displaying Only the Items of Interest in a Result Table. Printing a Report Using a Defined Template. Starting a New Study and Setting Up an Analysis About Studies and Analyses. Studies.	
Chapter 3	Databases Setting Up the Expected Compounds Library for a Targeted Analyses Checking the Access to the Online Databases for Identifying Unknown Compounds Setting Up a Study and an Analysis and Processing a Data Set Setting Up a New Study and a New Analysis by Using the Wizard Troubleshooting an Analysis Submitting a Run to the Job Queue Opening the Result File, Reviewing the Data, and Creating a Report Opening the Result File from the Job Queue Reviewing the Compounds that the Analysis Detected or Found. Displaying Only the Items of Interest in a Result Table. Printing a Report Using a Defined Template. Printing a Run Study and Setting Up an Analysis About Studies and Analyses. Studies. Input Files.	
Chapter 3	Databases Setting Up the Expected Compounds Library for a Targeted Analyses Checking the Access to the Online Databases for Identifying Unknown Compounds Setting Up a Study and an Analysis and Processing a Data Set Setting Up a Study and an Analysis and Processing a Data Set Setting Up a New Study and a New Analysis by Using the Wizard Troubleshooting an Analysis Submitting a Run to the Job Queue Opening the Result File, Reviewing the Data, and Creating a Report Opening the Result File from the Job Queue Reviewing the Compounds that the Analysis Detected or Found. Displaying Only the Items of Interest in a Result Table. Exporting a List of Compounds to a Spreadsheet Printing a Report Using a Defined Template. Starting a New Study and Setting Up an Analysis About Studies and Analyses. Studies. Input Files. Study Factors.	
Chapter 3	Databases Setting Up the Expected Compounds Library for a Targeted Analyses Checking the Access to the Online Databases for Identifying Unknown Compounds Setting Up a Study and an Analysis and Processing a Data Set Setting Up a New Study and a New Analysis by Using the Wizard Troubleshooting an Analysis Submitting a Run to the Job Queue Opening the Result File, Reviewing the Data, and Creating a Report Opening the Result File from the Job Queue Reviewing the Compounds that the Analysis Detected or Found. Displaying Only the Items of Interest in a Result Table. Printing a Report Using a Defined Template. Printing a Run Study and Setting Up an Analysis About Studies and Analyses. Studies. Input Files.	

	Data Storage Architecture	70
	Starting the New Study and Analysis Wizard	71
	Setting Up the Study Folders and Selecting a Processing Workflow	72
	Naming a Study and Its Folder	73
	Creating or Selecting the Top-Level Folder for the Study	73
	Selecting a Study Template	74
	Selecting a Processing Workflow for the Analysis	74
	Adding Input Files to a New Study	75
	Characterizing New Input Files	77
	Selecting the Delimiters for Parsing the File Names	78
	Adding Categorical Study Factors	78
	Adding Numeric Study Factors	80
	Adding Biological Replicate Study Factors	81
	Deleting Study Factors	82
	Duplicating Study Factors	82
	Editing Study Factors	83
	Automatically Assigning Study Factor Values	83
	Manually Selecting the Study Factor Values	
	Resetting the Sample Assignments	84
	Selecting the Sample Types	84
	Using the Regular Expression Builder to Extract the Study Factors	85
	Setting Up the Sample Groups and Ratios for a New Analysis	
	Opening the Wizard's Sample Groups and Ratios Page	92
	Selecting the Study Variables for Sample Grouping	
	Fixing Assignment Errors for Unset Study Variables	
	Changing the Hierarchy of the Study Variables	95
	Individually Setting Up the Group Ratios	97
	Using the Bulk Ratio Generation Feature to Set Up the Group Ratios .	97
	Preparing to Submit a Run to the Job Queue	
	About Accessing the Wizard Help	98
Chapter 4	Editing Existing Studies	99
•	About Existing Study Files	
	Opening an Existing Study	
	Study Page Commands and Tabs	
	Editing the Study Factors and the Study Description	
	Editing the Study Description	
	Adding, Deleting, or Editing Study Factors in an Existing Study	
	Study Definition Page of a Study	
	Adding Input Files to an Existing Study	
	Removing Input Files or Updating Their Location	
	Displaying the Location of a Study's Input Files	
	Removing Input Files from a Study	
	Resolving the Location of a Study's Input Files	
	Input Files Page Parameters	

	Editing the Sample Type and Study Factor Values	
	Editing the Sample Identifier on the Samples Page	
	Editing the Sample Types on the Samples Page	111
	Editing the Study Factor Values on the Samples Page	111
	Viewing the File Information for Study Samples on the Samples Page	111
	Samples Page Parameters	112
	Saving a Study File	113
Chapter 5	Creating, Running, and Reprocessing Analyses	115
	Setting Up a New Analysis from Within an Existing Study	115
	Starting a New Analysis from Within an Existing Study	
	Selecting the Input Files for a New Analysis	118
	Naming the Result File for a New Analysis	
	Setting Up the Sample Groups and Ratios for a New Analysis	
	Troubleshooting Common Analysis Errors	
	Common Analysis Errors.	
	Submitting an Analysis to the Job Queue	
	Common Validation Issues	
	Working with the Job Queue	
	Opening the Job Queue Page	
	Pausing a Run That Is Waiting to be Processed	
	Resuming a Paused Run	
	Canceling a Run	
	Removing Completed or Failed Runs from the Job Queue	
	Refreshing the Job Queue	
	Opening a Result File from the Job Queue	
	Displaying Verbose Messages in the Job Queue.	
	Viewing the Processing Steps for a Job.	
	Filtering the Job Queue.	
	Job Queue Page Parameters	
	Reviewing or Reprocessing an Analysis	
	Analysis Results Page Parameters	
	Reviewing an Analysis	
	Reprocessing an Analysis	
	Analysis Pane Parameters	
Chapter 6	Creating and Editing Processing Workflows	135
Shaptor V	Processing Workflows	
	Workflows Page.	
	Opening the Workflows Page	
	Workflows Page Command Bar	
	Workflows Page Shortcut Menu	138

	Editing a Processing Workflow	139
	Fixing a Workflow Node That Has a Caution Symbol	139
	Fixing a Workflow Node that Has an Exclamation Mark	140
	Adding a Workflow Node to a Processing Workflow	140
	Deleting a Workflow Node from a Processing Workflow	141
	Editing the Parameter Settings for a Workflow Node	141
	Setting Up Isotope Patterns for Pattern Scoring or Pattern Traces	142
	Opening the Isotope Ratio Editor	142
	Defining an Isotope Pattern from an Elemental Composition	144
	Copying an Elemental Composition from the Expected Compounds	
	Library	
	Setting Up a Custom Isotope Pattern	
	Exporting a Mass Spectrum from a RAW File to the Clipboard	
	Isotope Ratio Editor Parameters	
	Creating an Isotope Patterns List	
	Extracting Analog Traces	
	Extracting PDA Traces	
	Creating a Completely New Processing Workflow	
	Connecting the Workflow Nodes	
	Input and Output Nodes for the Workflow Nodes	
	Saving a Processing Workflow as a Template	160
a -		100
Chanter /	Workflow Nodes	10.5
Chapter 7	Workflow Nodes	
Chapter /	Input and Output Nodes	163
Chapter /	Input and Output Nodes	163 164
Chapter /	Input and Output Nodes	163 164 165
Chapter /	Input and Output Nodes. Export Spectra Node. Input Files Node. Spectrum Processing Nodes.	163 164 165 165
Chapter /	Input and Output Nodes.Export Spectra NodeInput Files NodeSpectrum Processing Nodes.Align Retention Times Node	163 164 165 165 166
Chapter /	Input and Output Nodes. Export Spectra Node. Input Files Node . Spectrum Processing Nodes. Align Retention Times Node. Filter By Mass Defect Node.	163 164 165 165 166 167
Chapter /	Input and Output Nodes.Export Spectra NodeInput Files NodeSpectrum Processing Nodes.Align Retention Times Node	163 164 165 165 166 167 170
Chapter /	Input and Output Nodes. Export Spectra Node . Input Files Node . Spectrum Processing Nodes. Align Retention Times Node . Filter By Mass Defect Node . Filter By Scan Event Node . Filter Centroids Node .	163 164 165 165 166 167 170 173
Chapter /	Input and Output Nodes. Export Spectra Node. Input Files Node . Spectrum Processing Nodes. Align Retention Times Node. Filter By Mass Defect Node. Filter By Scan Event Node .	163 164 165 165 167 167 170 173 174
Chapter /	Input and Output Nodes. Export Spectra Node. Input Files Node . Spectrum Processing Nodes. Align Retention Times Node. Filter By Mass Defect Node. Filter By Scan Event Node . Filter Centroids Node . Select Spectra Node.	163 164 165 165 166 167 170 173 174 181
Chapter /	Input and Output Nodes. Export Spectra Node . Input Files Node . Spectrum Processing Nodes. Align Retention Times Node. Filter By Mass Defect Node. Filter By Scan Event Node . Filter Centroids Node . Select Spectra Node . Trace Creation Nodes .	163 164 165 165 166 167 170 173 174 181 182
Chapter /	Input and Output Nodes.Export Spectra NodeInput Files NodeSpectrum Processing Nodes.Align Retention Times Node.Filter By Mass Defect Node.Filter By Scan Event NodeFilter Centroids NodeSelect Spectra Node.Trace Creation NodesCreate Analog Trace Node	163 164 165 165 166 167 170 173 174 181 182 184
Chapter /	Input and Output Nodes. Export Spectra Node. Input Files Node . Spectrum Processing Nodes. Align Retention Times Node. Filter By Mass Defect Node. Filter By Scan Event Node . Filter Centroids Node . Select Spectra Node. Trace Creation Nodes . Create Analog Trace Node .	163 164 165 165 166 167 170 173 174 181 182 184 186
Chapter /	Input and Output Nodes. Export Spectra Node . Input Files Node . Spectrum Processing Nodes. Align Retention Times Node. Filter By Mass Defect Node. Filter By Scan Event Node . Filter Centroids Node . Select Spectra Node . Trace Creation Nodes . Create Analog Trace Node . Create FISh Trace Node . Create Mass Trace Node .	163 164 165 165 166 167 170 173 174 181 182 184 186 187
Chapter /	Input and Output Nodes. Export Spectra Node. Input Files Node . Spectrum Processing Nodes. Align Retention Times Node. Filter By Mass Defect Node. Filter By Scan Event Node . Filter Centroids Node . Select Spectra Node . Create Analog Trace Node . Create FISh Trace Node . Create Mass Trace Node . Create Mass Trace Node .	163 164 165 165 166 167 170 173 174 181 182 184 186 187 189
Chapter /	Input and Output Nodes. Export Spectra Node . Input Files Node . Spectrum Processing Nodes. Align Retention Times Node. Filter By Mass Defect Node . Filter By Scan Event Node . Filter Centroids Node . Select Spectra Node . Create Analog Trace Node . Create FISh Trace Node . Create FISh Trace Node . Create Pattern Trace Node . Create Pattern Trace Node .	163 164 165 165 166 167 170 173 174 181 182 184 186 187 189 189
Chapter /	Input and Output Nodes. Export Spectra Node . Input Files Node . Spectrum Processing Nodes. Align Retention Times Node. Filter By Mass Defect Node. Filter By Scan Event Node . Filter Centroids Node . Select Spectra Node . Create Analog Trace Node . Create FISh Trace Node . Create FISh Trace Node . Create Pattern Trace Node . Create Pattern Trace Node . Compound Detection Nodes .	163 164 165 165 166 167 170 173 174 181 182 184 186 187 189 189 191
Chapter /	Input and Output Nodes. Export Spectra Node . Input Files Node . Spectrum Processing Nodes. Align Retention Times Node. Filter By Mass Defect Node. Filter By Scan Event Node . Filter Centroids Node . Select Spectra Node . Trace Creation Nodes . Create Analog Trace Node . Create FISh Trace Node . Create FISh Trace Node . Create Pattern Trace Node . Create Pattern Trace Node . Compound Detection Nodes . Analyze Labeled Compounds Node .	163 164 165 165 166 167 170 173 174 181 181 182 184 186 187 189 189 191 193
Chapter /	Input and Output Nodes. Export Spectra Node . Input Files Node . Spectrum Processing Nodes. Align Retention Times Node. Filter By Mass Defect Node . Filter By Scan Event Node . Filter Centroids Node . Select Spectra Node . Create Analog Trace Node . Create Analog Trace Node . Create FISh Trace Node . Create Mass Trace Node . Create Pattern Trace Node . Compound Detection Nodes . Analyze Labeled Compounds Node . Detect Compounds Node . Fill Gaps Node . Group Compounds Node . Peak Area Refinement Nodes .	163 164 165 165 166 167 170 170 173 174 181 182 184 186 187 189 189 191 193 195 196
Chapter /	Input and Output Nodes. Export Spectra Node . Input Files Node . Spectrum Processing Nodes. Align Retention Times Node. Filter By Mass Defect Node. Filter By Scan Event Node . Filter Centroids Node . Select Spectra Node . Create Analog Trace Node . Create Analog Trace Node . Create FISh Trace Node . Create Mass Trace Node . Create Pattern Trace Node . Create Pattern Trace Node . Compound Detection Nodes . Analyze Labeled Compounds Node . Fill Gaps Node . Group Compounds Node .	163 164 165 165 166 167 170 170 173 174 181 182 184 186 187 189 189 191 193 195 196

Compound Identification Nodes	200
Assign Compound Annotations Node	200
Predict Compositions Node	202
Search ChemSpider Node	206
Search Mass Lists Node	207
Search mzCloud Node	209
Search mzVault Node	214
Pathway Mapping Nodes	217
Map to BioCyc Pathways Node	217
Map to KEGG Pathways Node	219
Map to Metabolika Pathways Node	220
Compound Scoring Nodes	221
Apply Spectral Distance Node	222
Apply mzLogic Node	223
Calculate Mass Defect Node	224
Compound Class Scoring Node	225
Generate Molecular Networks Node (Beta)	226
Pattern Scoring Node	229
Expected Compounds Nodes	230
Find Expected Compounds Node	230
FISh Scoring Node	232
Generate Expected Compounds Node	234
Group Expected Compounds Node	239
Mark Background Compounds Node–Expected Compounds	240
Merge Features Node	241
Post-Processing Nodes	242
Descriptive Statistics Node	242
Differential Analysis Node	243
Export Xcalibur Inclusion or Exclusion List Node	243
Scripting Node	244
Reviewing the Analysis Results	247
Opening, Closing, and Updating Result Files	248
Opening Result Files Created in the Current Application Version	248
Opening Result Files Created in Previous Versions of the Application	249
Update Modes for Legacy Result Files	250
Closing a Result File	
Factory Default Layout for a Result Page	251
	Assign Compound Annotations Node

Modifying the Result Page Layout	253
Floating a Result Page View	253
Enlarging a Result Page View	253
Collapsible Pane Options for Filtering, Grouping, Coloring, and	
Discriminating By	253
Using a View's Collapsible Pane to Filter the Data By the Study	
Variable Values	
Using a View's Collapsible Pane to Filter the Data By Input File	256
Using a View's Collapsible Pane to Group the Data	256
Showing or Hiding Result Tables	257
Saving, Restoring, and Managing Layouts	257
Saving the Current Layout of a Result File	258
Resetting the Layout to the Factory Defaults	258
Creating a Custom Layout	258
Applying a Custom Layout	259
Managing the Layouts	259
Editing Compound Annotations	259
Adding and Deleting Proposed Structures for a Compound	261
Adding Structure Proposals	262
Deleting Structure Proposals	262
Replacing an Annotation with a Structure Proposal	262
Applying FISh Scoring	262
Applying FISh Scoring from a Result Table	263
Applying FISh Scoring from the Compound Annotation Editor Dialog	
Box	264
Specifying the FISh Scoring Parameter Settings	264
Using Result Filters for Data Reduction and Creating Filter Sets	266
Setting Up, Applying, and Saving Filter Sets	266
Creating a Result Filter for a Status Column	269
Creating a Result Filter with an AND Logical Conjunction	271
Creating a Result Filter with an OR Logical Conjunction	271
Creating a Result Filter with Both of the Logical Conjunctions	272
Loading a Saved Filter Set	273
Result Filters View Parameters	274
Viewing the Result Summary	275
Workflow Summary	276
Processing Messages Summary	276
Filter Summary	276
Study Summary	276
Grouping & Ratios Summary	276

	Exporting the Tabular Results to an External File	277
	Shortcut Menu Commands for the Result Tables	
	Exporting the Result Table Contents to a Spreadsheet	
	Exporting the Result Table Contents to a Text File	
	Exporting an Xcalibur Inclusion or Exclusion List	
	Exporting the Contents of the Compounds Table to TraceFinder	
	Exporting Spectral Data to a New or Existing mzVault Library	
	Adding a Compound to an Existing mzVault Library	
	Creating a New mzVault Library	
	Exporting Compounds to a New or Existing Mass List	
	Exporting Compounds from the Compounds Table to a New Mass	
	List	289
	Exporting Compounds from the Compounds Table to an Existing	
	Mass List	
	Copying or Saving Graphical Views for Publication	291
Chapter 9	Working with the Graphical Result Views	
	Working with the Chromatograms View	
	Viewing a Chromatogram	294
	Adding a Chromatogram Plot	
	Overlaying Chromatogram Plots	
	Changing the Grouping in the Collapsible Pane	
	Hiding the Traces for a Study Variable Value	
	Updating All the Chromatogram Plots Simultaneously	
	Manually Integrating Chromatographic Peaks	
	Chromatograms View Shortcut Menu Commands	
	Working with the Mass Spectrum View	305
	Displaying a Mass Spectrum	
	Changing the Zoom Level of the Mass Spectrum View	
	Viewing Fragment Structures for Expected Compounds	
	Creating a Mirror Plot	307
	Searching the mzCloud Library for a Matching Fragmentation	
	Spectrum	
	Spectral Tree Pane of the Mass Spectrum View	
	Isotope Pattern Matching for Compounds with Formulas	
	Mass Spectrum View Shortcut Menu Commands	
	Working with the Result Charts	
	Opening the Result Charts View	312
	Using the Copy, Export, and Zoom Commands for a Result Charts View.	312
	Displaying and Pinning the Options Pane for the Result Charts View	
	Plotting the Data as a Histogram Chart.	
	Plotting the Data as a Bar Chart	
	Plotting the Data as a Pie Chart.	
	I IOILIIIZ UIC D'ALA AS A I IC OIIAIL	

Working with a Scatter Plot	. 326
Setting Up a Scatter Plot	. 327
Using a Filter Set to Filter the Scatter Plot	. 328
Customizing the Appearance of a Scatter Plot	. 328
Customization Options for a Scatter Plot	. 329
Scatter Chart Parameters	. 330
Scatter Chart Shortcut Menu Commands	. 330
Working with the Trend Chart View	. 332
Opening the Trend Chart View	. 334
Defining the Sample Groups to Compare	. 334
Comparing the Peak Areas for a Single Compound by Sample Group	. 334
Changing the Sort Order of the Defined Groups	. 335
Comparing the Peak Areas for Multiple Compounds by Group	. 336
Changing the Hierarchy of the Variables Used for Grouping	. 337
Showing the Error Bars in a Trendline Chart	. 337
Trend Chart View Parameters	. 339
Working with the Mass Defect Plot View	. 339
Working with the Principal Component Analysis View	. 341
Setting Up a Principal Component Analysis	. 341
Interpreting the Scores Plot	. 342
Interpreting the Loadings Plot	. 343
Working Interactively with the Loadings Plot	. 344
Interpreting the Variance Plot	. 345
Principal Component Analysis Parameters	. 347
Working with the Descriptive Statistics View	. 347
Working with the Differential Analysis View.	. 352
Reviewing the Initial Differential Analysis	
Changing the Analysis Settings for a Differential Analysis	. 353
Running a New Differential Analysis	. 354
Differential Analysis View Parameters and Shortcut Menu Commands	. 356
Working with the Partial Least Squares Discriminant Analysis View	
Identifying a Set of Compounds to Discriminate Between Groups	
Partial Least Squares - Discriminant Analysis View Parameters	. 359
Viewing KEGG Pathways	. 360
Viewing BioCyc Pathways	. 362
Viewing Metabolika Pathways	
Viewing the Corrected Retention Times of the Alignment Features	. 365
Viewing the QC-Based Compound Area Correction Curves	. 367
Viewing the FISh Scoring Queue	. 368
Running a Hierarchical Clustering Analysis	. 369
Setting Up a Hierarchical Clustering Analysis	
Working Interactively with the Heat Map Grid	
Hierarchical Cluster Analysis View Parameters	. 373

	Running an mzLogic Analysis	
	Performing an mzLogic Analysis	
	Reviewing the Results of an mzLogic Analysis	
	mzLogic Analysis View Parameters	
	Working with the Isotopologues Distribution Chart	379
Chapter 10	Reviewing the Tabular Result Data	
	Common Result Tables	
	Adducts Table	
	Chromatogram Peaks Table	
	File Alignments Table	
	FISh Trace Fragments Table	
	Input Files Table	
	Manual Peaks Table	
	Merged Features Table	
	Specialized Traces Table	
	Structure Proposals Table	
	Expected Compounds Result Tables	
	Expected Compounds Table	
	Expected Compounds per File Table	
	Expected Features Table	
	Expected Formulas Table	
	Related Structures Table	
	Transformations Table	
	Compound Detection Result Tables	
	Compounds Table	
	Compounds per File Table	
	Features Table	
	Filled Gaps Table	
	Labeled Features Table	
	Labeled Compounds per File Table	
	Similar Compounds (Beta) Related Table	
	Compound Identification Result Tables	
	ChemSpider Results Table	
	Mass List Search Results Table	
	mzCloud Results Table	
	mzVault Results Table	
	Predicted Compositions Table	
	Pathway Mapping Result Tables	
	BioCyc Pathways Table	425
	BioCyc Results Table	
	KEGG Pathways Table	
	Metabolika Pathways Table	
	Metabolika Results Table	

	Compound Scoring Tables
	Compound Class Matches Table
	Matched Patterns Table
	Differential Analysis Columns
	Descriptive Statistics Columns
Chapter 11	Creating and Printing Reports437
	Reporting Workflow
	Generating a Report with an Existing Report Template
	Creating a New Report Template
	Creating a New Report Template by Using the Customize Report
	Dialog Box
	Adding, Removing, or Modifying Color Schemes for a Report
	Template's Tables
	Customize Report Dialog Box Parameters
	Editing an Existing Report Template
	Opening a Report Template
	Changing a Report Template's Logo Image
	Changing a Report Template's Time Stamp
	Adding Design Items to a Report Template
	Adding Main Table Columns to a Report Template
	Adding Data Graphs to a Report Template
	Adding Related Table Columns to a Report Template
	Editing the Properties of Subreport Columns in a Report Template458
	Moving a Subreport Column to a Report Template's Main Table459
	Modifying the Properties of a Report Template Design Item
	Adding Page Breaks to a Report Template
	Deleting a Pair of Workspace Sections on the Report Template Page 462
	Resizing a Report Template's Sections
	Editing Tools for the Report Template Page
	Workspace Sections and Sizing Bar on the Report Template Page
	Report Template Page Toolbars
	Report Template Page Shortcut Menu
	Section Report Items for a Report Template
	Property Settings for the Sections and Items in a Report Template
	Appearance Properties
	Behavior Properties
	Data Properties
	Design Properties
	Layout Properties
	Miscellaneous Properties
	Opening the Property Dialog Boxes for the Report Template's Design
	Items
	Populating a Rich Text Box by Loading a File

	Selecting the Paper Type, Print Width, Page Orientation, and Watermark	
	for a Report Template	487
	Previewing and Printing a Report	489
	Report Preview Toolbar	493
	Page Thumbnails Pane	495
Chapter 12	Using the Lists and Libraries Manager	497
	Library Manager Page	497
	Modifying the Expected Compounds List	498
	Importing, Exporting, and Deleting Compounds	498
	Adding and Editing Compounds with the Compound Editor	500
	Modifying the Adducts List	502
	Working with the Adducts List	503
	Adding and Editing Adducts with the Adduct Editor	503
	Modifying the Ion Definitions List	505
	Working with the Ion Definitions List	
	Adding or Editing Ion Definitions with the Ion Definition Editor	508
	Modifying the Transformations List	510
	Deleting, Importing, or Exporting Transformations	510
	Adding or Editing Transformations with the Transformation Editor	
	Modifying the Mass Lists Library	
	Adding, Deleting, Importing, Exporting, and Replacing Mass Lists	518
	Editing a Mass List	521
	Adding and Editing Mass List Compounds with the Compound Editor .	
	Modifying the Spectral Libraries List	
	Modifying the Metabolika Pathways List.	
	Deleting, Importing, Exporting, and Replacing Metabolika Pathways	
	Editing Metabolika Pathways	
	Adding and Editing Pathway Structures with the Compound Editor	
	Modifying the Compound Classes List	
	Compound Class Editor	
	Using the Fragment Editor	
	Loading a Structure from a Structure File	
	Finding a Structure in the ChemSpider Database	
	Using the Structure Drawing Tools or Commands	
	Compound Editor Toolbar and Fragment Editor Toolbar	
	Shortcut Menu Commands for the Drawing Area	
	Using the Structure Icons	
	Using the Template Tool	
	Checking the Validity of a Structure	
	Manipulating Structures	
	Modifying Atoms and Bonds.	
	Saving a Structure	
	Editing Atom Properties	550

Chapter 13	Using the License Manager	
	Opening the License Manager	553
	License Manager Command Bar	554
	Activating the Software License	555
	Entering the Product ID and the Activation Code	555
	Activating the License on an Online Computer	557
	Activating the License on an Offline Computer.	
	Deactivating the Software License for Transfer to Another Computer	560
	Deactivating the Software License on an Online Computer	
	Deactivating the Software License on an Offline Computer	
	Installing or Updating a Processing Workflow Node	563
Chapter 14	Setting Up the Configuration Options	565
	Opening the Configuration Page	565
	Selecting the Maximum Number of Parallel Processing Jobs	565
	Selecting Where to Store Temporary Data	566
	Turning Off the Auto-Save Feature for Studies	567
	Hiding the Workflow Node Numbers	567
	Selecting the Global Color Palettes	568
	Opening the Color Schema View	569
	Selecting a Standard Color Palette	569
	Creating Custom Color Palettes	569
	Deleting Custom Color Palettes	570
	Importing Custom Color Palettes	570
	Exporting Custom Color Palettes	571
	Editing a Custom Color Palette	571
	Selecting a Color in the Gradient Color Chart	573
	Specifying the Default mzCloud Mass Tolerance Settings	574
	Setting Up a BioCyc Account or Subscription	575
	Opening the BioCyc User Login View	575
	Setting Up a BioCyc Account or Individual Subscription	576
	Entering, Testing, and Saving Your Account Information	576
	Specifying the Fragmentation Libraries	577
.		
Chapter 15	Working with the Application Tables	
	Moving Up and Down Table Rows.	
	Sorting Data Tables.	
	Sorting Table Entries By One or More Columns.	
	Sorting Table Entries By a Column with a Distribution Map	
	Freezing Table Rows	
	Grouping Table Rows	
	Changing the Position of Table Columns	
	Changing the Column Order	
	Stacking Two Table Columns into One Column	
	Freezing Table Columns	585

	Showing or Hiding Table Columns	585
	Copying Table Entries to the Clipboard	586
	Copying the Contents of a Single Table Cell to the Clipboard	586
	Copying the Contents of a Single Row to the Clipboard	
	Copying the Contents of Multiple Rows to the Clipboard	
	Filtering the Tables on a Study Page or a List or Library View	
	Setting Up Single-Condition Filters for the Table Columns	
	Setting Up a Single-Condition Wild Card Filter	
	Setting Up a Single-Condition Filter for Numeric Data	589
	Operators and Operands for a Single-Condition Table Filter	589
	Setting Up a Custom Filter with Multiple Conditions	594
	Custom Filter Selection Dialog Box	598
Chapter 16	Working with Molecular Networking	
•	Using the Molecular Networking Feature	
	How the Generate Molecular Networks Node Works	600
	Information Displayed in the Similar Compounds Table	
	Exporting the Molecular Network to the Viewer	
	Working with the Molecular Network Viewer.	
	Color-Coding for the Nodes	
	Sizing a Node By the Compound's Peak Area or MW	
	Colorizing a Link By Its Score, Coverage, or Number of Fragments .	605
	Interactive Functions in the Molecular Network Viewer	606
	Search Pane	606
	Graph Info Pane	607
	Isolation Pane	607
	Filters Pane	607
	Thresholds Pane	608
	Compound or Transformation Information	
	Using the Isolation mode	609
Chapter 17	Testing Communication to the Online Databases	611
-	Running the Communication Tests	
	Checking the URLs for the Online Databases in Your Browser	
	Specifying the IP Address of the Proxy Server	614
	Setting the Correct Time and Time Zone on the Processing Computer	615
Appendix A	Experiment Design for Comparison Statistics	617
	Biological Versus Technical Replicates	
	Non-Nested Versus Nested Experiment Designs	
	Index	

Preface

This guide describes how to use the Compound Discoverer[™] application to qualitatively process RAW data files with a targeted or untargeted workflow. A targeted workflow evaluates the mass spectral data for the presence of specific compounds. An untargeted workflow evaluates the mass spectral data, predicts the elemental composition of unknown compounds, and searches mass spectral databases to identify these compounds.

To familiarize yourself with the Compound Discoverer application, follow the tutorials that are available from the application Help menu or see Chapter 2, "Getting Started."

Contents

- Related Documentation
- System Requirements
- Installation Instructions
- Special Notices
- Contacting Us

Related Documentation

The Compound Discoverer application includes these manuals as PDF files:

- Compound Discoverer User Guide
- Compound Discoverer E & L Tutorial
- Compound Discoverer Metabolism Tutorial
- Compound Discoverer Metabolomics Tutorial
- Compound Discoverer Stable Isotope Labeling Tutorial
- Compound Discoverer Reporting Quick Start

The Compound Discoverer application also includes a Help system.

* To view the Compound Discoverer manuals

From the application window, choose **Help > Manuals**.

-or-

From the Microsoft[™] Windows[™] taskbar, choose **Start > All Programs (or Programs) > Thermo Compound Discoverer 3.1 >** *Manual*.

★ To download user documentation from the Thermo Scientific[™] website

- 1. Go to thermofisher.com.
- 2. Point to Support, and then click Manuals under Product Documentation on the left.
- 3. In the Refine Your Search box, search by the product name.
- 4. From the results list, click the title to open the document in your web browser, save it, or print it.

To return to the document list, click the browser **Back** button.

System Requirements

The Compound Discoverer 3.1 application can process data files produced by high-resolution accurate-mass (HRAM) Thermo Scientific[™] mass spectrometers, such as the Orbitrap Fusion[™], Q Exactive[™], and Exactive[™].

Table 1 lists the hardware and software requirements for the processing computer.

Table 1. Hardware and software requirements for the processing computer

System	Minimum requirements
Hardware	• 3.4 GHz dual-core processor
	• 16 GB RAM
	• 500 GB hard drive
	• DVD-ROM and USB drive
	• Display monitor resolution of 1920 × 1080 with 96 dpi setting
Software	 Microsoft Windows 7 Pro SP1 (64-bit) or Windows 10 64-bit operating system
	• Microsoft .NET Framework 4.7.2
	Microsoft Office 2010
	 Adobe[™] Reader[™] 11
	• Adobe Flash™ Player 15
System settings	• To run processing workflows with online mass spectral database searches, the computer must have unblocked access to the mass spectral databases on the Internet.
	• The computer must have the correct time and date settings and be synchronized with Internet time.
	• The Region and Language setting for the operating system must be set to English (United States).

Table 2 lists the recommended hardware configurations for enhanced performance using theCompound Discoverer application.

System	Recommended configurations
Hardware	 Dual 8-core processor (for example, 2x Intel[™] Xeon[™] Gold 6134 CPU @ 3.20 GHz)
	• 64 GB RAM
	• 1 TB SSD (solid-state disk) hard drive for OS
	• 2nd 3 TB (conventional disk) hard drive for data storage
	• DVD-ROM and USB drive
	 Two 27 in. UHD monitors: Display monitor resolution of 3840 × 2160

Table 2. Recommended hardware configurations for enhanced performance

To check the access to the mass spectral databases, the time and date settings, and the Internet time, see Chapter 17, "Testing Communication to the Online Databases."

To verify that the system meets the minimum requirements, follow these procedures:

- To check the computer specifications
- To check the font DPI for a Windows 7 system
- To check the format setting for Region and Language

To check the computer specifications

1. From the Windows Explorer directory, right-click **OSDisk** (*Drive*:) (the directory for the hard drive where the operating system is installed) and choose **Properties**.

The OSDisk (Drive:) Properties dialog box opens. This dialog box lists the file system (NTFS or FAT) and the free disk drive space.

From the Windows Desktop, choose Start > Control Panel > System and Security > System.

The System page opens. This page lists the operating system; the processor type, speed, and number of cores; the installed RAM; and the system type (32-bit or 64-bit).

To check the font DPI for a Windows 7 system

- 1. From the Windows taskbar, choose **Start > Control Panel**.
- 2. On the Adjust Your Computer's Settings page, do the following:
 - a. For View By, select **Category**.

- b. Under Appearance and Personalization, click Adjust Screen Resolution.
 The Screen Resolution page opens.
- 3. Click Make Text and Other Items Larger or Smaller.
- 4. On the left panel, click Set Custom Text Size (DPI).

The Custom DPI Setting dialog box opens (Figure 1).

Figure 1. Custom DPI Setting dialog box

Custom DPI Sett	ting		x
	OPI setting, select a with your mouse.	percentage from t	he list, or
Scale to this pe	ercentage of norma	size: 100% 🔻	
0	1	2	3
9 point Sego	e UI at 96 pixels pe	r inch.	
√ Use Windo	ws <u>X</u> P style DPI scal	ing OK	Cancel

5. Make sure that the DPI setting is 96 pixels per inch.

To check the format setting for Region and Language

- 1. From the Windows taskbar, choose **Start > Control Panel**.
- 2. On the Adjust Your Computer's Settings page, do the following:
 - a. For View By, select **Category**.
 - b. Select Clock, Language, and Region.
- 3. Select Region and Language.
- 4. On the Formats page of the Region and Language dialog box, select **English** (**United States**) from the Format list.

Installation Instructions

Thermo Compound Discoverer is a licensed application. After you install the application, you can use it for up to 60 days without activating the software license.

After you order the Compound Discoverer 3.1 application, you will receive a software media kit that includes a key-shaped USB flash drive with the installation executable. In addition, you will receive an email from Thermo MS Licensing providing you with the information that you need to activate the software license.

If you are upgrading the software from a previous version of the application, you can find the software installer and a license on the LSMS Software Download and Licensing Portal.

IMPORTANT Read the following:

- The installation requires the Windows 7 SP1 64-bit or Windows 10 64-bit operating system. See "System Requirements" on page xix for the recommended hardware requirements and system settings.
- The Compound Discoverer 3.1 licensing process requires an Internet connection to validate the software license. You can install the application on a computer without Internet access and complete the activation process on a computer with Internet access.
- The Compound Discoverer application is supported for US-English Only locale.

Note The following versions of the Compound Discoverer application can coexist on the same computer: 1.0, 2.0, 2.1, 3.0, and 3.1

The installation executable includes three installers. The Thermo Compound Discoverer 3.1 installer installs the Compound Discoverer application. The Thermo mzVault Library installer installs two mzVault libraries that are May 2019 snapshots of the online mzCloud mass spectral database, and the Thermo mzVault installer installs the mzVault application.

To install the applications and the libraries, see these topics:

- Installing the Compound Discoverer Application and the mzVault Libraries
- Installing the mzVault 2.2 Application

Installing the Compound Discoverer Application and the mzVault Libraries

* To install the Compound Discoverer application and the mzVault libraries

- 1. Do one of the following:
 - a. Insert the Compound Discoverer USB flash drive into a USB port on your computer.
 - b. Open Windows Explorer and select the USB drive to view its contents.

c. Locate the executable: XStart_Compound Discoverer.exe.

-or-

a. Go to the following website address:

https://thermo.flexnetoperations.com

- b. Log in to your account. If you do not have an account, click **Register** and create one.
- c. On the left side of the Life Sciences Mass Spectrometry Software Download and Licensing Portal, click **Product List** under Software and Services.
- d. On the Product List page, click the **Demo** hyperlink.
- e. On the Product Information page, click the **Compound Discoverer 3.1** hyperlink.
- f. On the Product Download page, in the File Name column, click the down arrow to the left of Compound Discoverer 3.1.zip, click **Save As**, and then save the compressed zipped folder to your computer.
- g. Extract the contents of the zipped folder, and locate the executable: XStart_Compound Discoverer.exe.

Note If you are upgrading from the Compound Discoverer 3.0 software, which is a free upgrade, you can also find your software license on the Product Download web page.

2. Double-click XStart_Compound Discoverer.exe.

The installation wizard starts.



Adobe Acrobat Reader or an alternative PDF reader must be installed before you can Adobe® Reader® XI Copyright © 1984-2012. Adobe® Flash® Player. Copyright © 19	read any of the documents provided as F 96-2010. Adobe Systems Incorporated. A	DF files on this media. Validation certificate is also included. Il Rights Reserved. Adobe Reader and Flash Player are eithe
trademarks or registered trademarks in the Unites States and/or other countries.		
	DMPOUND DISCOVERER	Installation Guide
	0 18-	Release Notes
		Compound Discoverer 3.1
		mzVault 2.2
Compound Discoverer 2	-	mzVault 2.2 Release Notes
Compound Discoverer 3. Integrated Solutions for Small Molecule Research		mzVault Library Installer
		Microsoft .NET Framework 4.7.2
2.02	4	Adobe Reader 11.0
© Copyright 2014-2019 Thermo Fisher Scientific Inc. All rights reserved. This program is protected by copyright law and international treaties as described in Help About.	scientific	Adobe Flash Player 15.0
		Browse Media (Example Studies,)

3. Click Compound Discoverer 3.1.

4. When the installer opens, click **Next** and follow the instructions.

The Minimum Product Requirements screen appears with a product requirements scan summary. If the computer meets the system requirements, the summary states the following:

All the system minimum requirements checks passed.

- 5. At the Installation Qualification prompt, click **Yes** if you want to view the Installation Qualification report. Otherwise, click **No**.
- 6. When the installation is complete, select whether you want the installer to restart your computer, and then click **Finish**.
- 7. To install the mzVault libraries, do the following:
 - a. If the installer restarted your computer, double-click **XStart_Compound Discoverer.exe** to restart the installer.
 - b. Click mzVault Library Installer to install the mzVault libraries:
- 8. If you have not already restarted your computer, restart it now.

Installing the mzVault 2.2 Application

To create mass spectral libraries from your own data acquired with a high-resolution accurate-mass (HRAM) Thermo Scientific mass spectrometer, install the mzVault 2.2 application. The Compound Discover installer includes an installer for the mzVault application.

To install the mzVault application

- 1. If the Compound Discoverer installer is not open, open it as follows:
 - a. Locate the XStart_Compound Discoverer executable.

You can find the executable on the USB drive provided in the Compound Discoverer application media or in the zip folder that you downloaded from the product download web site.

- b. Double-click XStart_Compound Discoverer.exe.
- 2. Click mzVault 2.2.

The Thermo mzVault Installshield Wizard opens.

Note If you have an earlier version of the application, follow the instructions to remove it, and then restart the mzVault 2.2 installer.

If your computer meets the minimum requirements, the Next button becomes available.

3. Click **Next** to continue.

- 4. When the installation is complete, do the following:
 - a. To start the mzVault application immediately after you exit the installer, select the **Launch Thermo mzVault** check box.
 - b. Click Finish.
 - c. If you want to view the Installation Qualification report, click **Yes** at the confirmation prompt. Otherwise, click **No**.

Special Notices

Make sure you follow the precautionary statements presented in this guide. The special notices appear in boxes.

Special notices include the following:

IMPORTANT Highlights information necessary to prevent damage to software, loss of data, or invalid test results; or might contain information that is critical for optimal performance of the system.

Note Highlights information of general interest.

Tip Highlights helpful information that can make a task easier.

Contacting Us

Contact	Email	Telephone	QR Code
U.S. Technical Support	us.techsupport.analyze@thermofisher.com	(U.S.) 1 (800) 532-4752	
U.S. Customer Service and Sales	us.customer-support.analyze@thermofisher.com	(U.S.) 1 (800) 532-4752	
Global support	 To find global contact information or custor 	nize your request	
	1. Go to thermofisher.com.		
	2. Click Contact Us , select the country, and the you need.	en select the type of support	
	3. At the prompt, type the product name.		
	4. Use the phone number or complete the onlin	e form.	
	 To find product support, knowledge bases, 	and resources	
	Go to thermofisher.com/us/en/home/technic	al-resources.	
	 To find product information 		
	Go to thermofisher.com/us/en/home/brands/	thermo-scientific.	
	back for this document, go to surveymonkey.com/s s (techpubs-lcms@thermofisher.com).	/PQM6P62 or send an email	message to

* For Compound Discoverer customer support questions

Send an email message to CD.support@thermofisher.com.

1

Introduction

Compound Discoverer is a qualitative data-processing application that uses accurate mass data, isotope pattern matching, and mass spectral library searches for the structural identification of small molecules. It can process the accurate-mass spectra from the entire product line of Thermo Scientific high-resolution mass spectrometers. It can also display the graphical data acquired from a variety of detectors: UV-visible and photodiode array (PDA) detectors that are controlled by a Thermo Scientific data system and third-party analog detectors that are connected to the analog input channels of a Thermo Scientific MS.

To familiarize yourself with the Compound Discoverer application, see these topics.

Contents

- New Features and Enhancements
- Supported File Formats
- Starting the Application
- Choosing the Toolbar Icon Size
- The Application Window
- Understanding Processing Workflows
- Best Scans for Composition Prediction and Spectral Matching
- Quality Control Samples for Batch Normalization
- Stable Isotope Labeling
- FISh Scoring for Targeted Compounds and Proposed Structures
- Calculating the Mass Defect of an Elemental Composition
- Using mzLogic to Score Candidates for Unknown Compounds
- Managing the Start Page's Recent File Lists
- Auto-Hiding the Start Page, the Chromatograms View, and the Mass Spectrum View
- Working with the Tabbed Documents
- Rearranging the Tabbed Documents and Graphical Views

Tip To get started with this application, follow these tutorials and guides or go to Chapter 2, "Getting Started."

- Compound Discoverer 3.1 Metabolism Tutorial
- Compound Discoverer 3.1 Metabolomics Tutorial
- Compound Discoverer 3.1 Extractables and Leachables Tutorial
- Compound Discoverer 3.1 Stable Isotope Labeling Tutorial
- Compound Discoverer 3.1 Reporting Quick Start Guide

To access these tutorials from the application window, choose Help > Manuals.

New Features and Enhancements

The application uses a study format to define the sample types, experimental study factors, sample groups, and group ratios. It uses a customizable node-based processing workflow to process Xcalibur[™] RAW files and create a result file.

Some of the workflow nodes require input from the application's customizable lists and libraries, which include the structures of known compounds, mass lists, adduct ions, transformations, metabolic pathways, fragment lists, and spectral libraries.

The result file includes a set of result tables and graphical views based on the processing workflow. The application provides graphical views for displaying chromatograms and spectra, and statistical views for comparing the detected compounds by sample file or sample group.

These topics describe the new features and enhancements:

- New Features
- Enhancements

New Features

Compound Discoverer 3.1 has the following new features:

- Molecular Networks (beta) to interactively explore relationships between compounds based on transformations and spectral similarity
- Scripting node to embed scripts in workflows—for example, Python or R scripts

Tip For information about using the Scripting Node, go to the Resources page of the following web site: https://mycompounddiscoverer.com/.

Mass Defect Plot for Compounds and Expected Compounds

- Calculate Mass Defect node with support for up to 5 Kendrick formulas
- Peak area scaling using values from a numerical study factor that you specify as the scaling factor in the Normalize Areas node
- Search mzCloud node:
 - Supports automatic searches of multiple spectral libraries
 - Supports the new "Autoprocessed" mzCloud library
 - Supports using both DIA and DDA scans as the query scans to search against the mzCloud library
 - Highlights matching and non-matching spectral peaks in green and red, respectively
 - Displays the fragment structure for matching peaks
- Apply Spectral Distance node to score the observed isotopic pattern versus the simulated isotopic pattern of a candidate from, for example, ChemSpider
- Manual export of detected compounds in a result file to a mass list file
- Installation of the mzVault 2.2 application for creating and editing mass spectral libraries

Note The mzVault 2.2 application supports importing NIST MSP and MassBank MB files.

Enhancements

Compound Discoverer 3.1 has the following enhancements:

- Improved performance of ChemSpider searches
- Displays the chromatographic peak areas in units based on a time scale of seconds instead of minutes
- Extended retention time search window for the assignment of MSn spectra to a compound
- Changes in the Expected Compounds database are now immediately reflected in the Generate Expected Compounds node
- New parameter for Compound Class Scoring node to use entire MSn tree
- New version of the HighChem Fragmentation Library
- New licensing mechanism

Supported File Formats

Table 3 describes the file types that the application can recognize or create.

 Table 3.
 Supported file formats (Sheet 1 of 2)

File format	Description
Xcalibur RAW file	Contains unprocessed data acquired from a high-resolution, accurate mass LC/MS/MS instrument with a Thermo Scientific data system that is layered on the Thermo Foundation [™] platform.
MOL format (.mol), compressed structure (.mcs), template (.tml)	Contains a two-dimensional compound structure. You can open structure files by using the Structure Editor or the Custom Explanations Editor.
cdProcessingWF	Contains the data processing instructions for the application. To create a processing workflow, you must start or open an analysis in a study.
cdAnalysis	Stores the processing workflow information.
cdStudy	Stores the study information, which includes the names and locations of the input files, the sample information, and the relationship between the input files.
cdResult	Contains the results produced by processing a set of raw data files and information about the analysis settings used to process the raw data.
cdResultView	Contains the layout settings that the application uses to display the result file's tables and graphical views. These settings also include the applied result filters. Deleting this file erases all the custom layout settings and restores the display to the default layout settings.
cdReportTemplate	Contains the layout for reports that extract data from the following items in a result file: selected columns in one main table, selected columns in the related tables, and selected graphical views (MS1, MS2, and Chromatograms).

File format	Description
Filter Set (.filterset)	You can use filter set files for data reduction when reviewing and reporting the data in Compound Discoverer result files. The application comes with one predefined filter set file—Example Filter Set.filterset.
mgf, mzML, mzDATA	Contains the mass spectral data that can be read by third-party mass spectrometry applications. To create any of these files types, use the Export Spectra Node in a processing workflow.
XML	You can export each library to an XML file, and you can import library entries from an XML file.
text (.txt)	You can save the data points in the graphical views to a text file.
EMF, BMP, JPG, GIF, PNG, TIFF	You can save the images in the graphical views as image files of the following file types: enhanced metafile (.emf), bitmap (.bmp), Joint Photographic Group (.jpg), graphic interchange format (.gif), portable network graphics (.png), and tagged image file format (.tiff). You can open EMF files in a raster image editor or a vector image
	editor.
CSV	You can import the contents of a CSV file into a Mass List file.
CLIB	You can import the contents of a CLIB file as a Compound Class list.
DB	You can import mzVault libraries into the Spectral Libraries list.

Table 3. Supported file formats (Sheet 2 of 2)

Starting the Application

You can start the application from the taskbar or the computer desktop.

- To start the application
 - From the taskbar, choose **Start > All Programs** (or **Programs**) **> Thermo Compound Discoverer**.

-or-

• From the computer desktop, double-click the **Compound Discoverer** icon,

The Compound Discoverer window opens with the Start Page displayed as a tabbed document (Figure 3). As you create studies and process data, the application creates and populates recent file lists to the right of the What Would You Like to Do? hyperlinks.

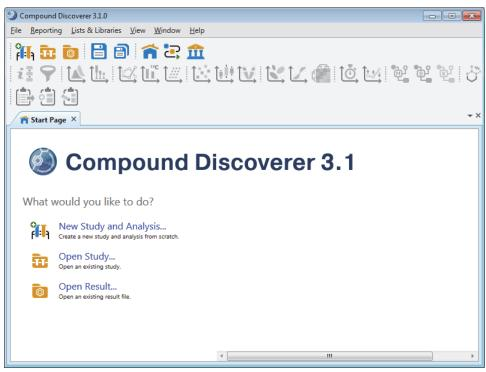


Figure 3. Application window with the initial Start Page and large toolbar icons

Choosing the Toolbar Icon Size

You can choose between large or small icons for the application toolbar. By default, the application toolbar displays large icons.

✤ To change the size of the toolbar icons

Do one of the following:

- To display small icons in the toolbar, right-click the application toolbar and choose ✓ Show Large Icons.
- To display large icons in the toolbar, right-click the application toolbar and choose **Show Large Icons**.

Note The check mark to the left of Show Large Icons indicates that the application is set to display large toolbar icons.

The Application Window

The application window contains a title bar, a menu bar, and a toolbar. From the application window, you can open all the other application views and pages by choosing a menu command or by clicking a toolbar icon.

Note This user guide uses the following terms to describe the user interface:

- View—A dockable window that you can move to a second monitor.
- Page—A tabbed document. You can have many pages open simultaneously; however, only one of these pages is the active page.
- Dialog box—A graphical element that accepts user input. Only one dialog box can be open at a time. When it is open, a dialog box blocks you from working in other parts of the application.
- Pane—A defined area of an application view, page, or dialog box.
- Prompt—A pop-up message box that you must dismiss to continue.

These topics describe the menu bar and the toolbar:

- Application Menu Bar
- Application Toolbar

Application Menu Bar

Table 4 describes the menu commands in the menu bar at the top of the application window.

File menu These commands are always available. New Study and Opens the New Study and Analysis Wizard, which takes you Analysis through the process of selecting the studies folder for your study subfolders, creating a new study, and starting a new analysis. Open Study Opens the Open Study dialog box for selecting an existing study file to open. The CD study file type has the .cdStudy file name extension.

Menu command	Description
Open Result	Opens the Open Result File dialog box for selecting an existing result file or result view file.
	You do not need an active software license to open result files.
	The CD result file type, which contains the data processing results, has the .cdResult file name extension.
	The CD result view file type, which contains the display layout for the results tables and graphical views and the filter settings, has the .cdResultView file name extension.
	To restore the default layout for a result file, delete its associated CD result view file.
	Note You do not need an active software license to open result files.
Save	Saves recent changes to the current active page (selected tab), for example, the active study page or result page.
Save All	Saves recent changes to all the open pages in the application window.
Recent Studies	Displays a list of recent studies that you can open.
Recent Studies > Clear	Clears the Recent Studies list.
Recent Results	Displays a list of recent results that you can open.
Recent Results > Clear	Clears the Recent Results list.

Table 4. Application menu bar (Sheet 2 of 6)

Reporting menu

These commands are only available when a result page is active.

Create Report	Opens the Open Report Design Template dialog box for selecting a report template to resolve specific data in the result file.
Create Report Template	Opens the Customize Report dialog box for setting up the main properties of a report template.
Edit Report Template	Opens the Open Report Design Template dialog box for choosing an existing report template to edit.

Table 4. Application menu bar (Sheet 3 of 6)

Menu command	Description	
List & Libraries menu		

These commands are always available.

Expected Compounds	Opens the Expected Compounds view for modifying the list of expected compounds.	
	Note The Generate Expected node and Create Pattern Trace node require compounds from the Expected Compounds library. The Create Pattern Trace and the Pattern Scoring nodes require user-specified elemental compositions.	
Transformations	Opens the Transformations view for modifying the list of transformations.	
Ion Definitions	Opens the Ion Definitions view for modifying the list of ion definitions.	
Adducts	Opens the Adducts view for modifying the list of adducts.	
Mass Lists	Opens the Mass Lists view for modifying the list of mass list files or editing mass list files.	
Spectral Libraries	Opens the Spectral Libraries view for modifying the list of mzVault database files.	
Metabolika Pathways	Opens the Metabolika Pathways view for modifying the list of Metabolika pathways or editing a Metabolika pathway.	
Compound Classes	Opens the Compound Class view for viewing or modifying the list of compound class libraries.	

View menu

The Start Page and Job Queue commands from this menu are always available. The other View commands are only available when a result file is active.

Start Page	Opens the Start Page, which lists the most recently opened result files and study files.
Job Queue	Opens the Job Queue page for viewing the progress of the current analysis or the processing events of previous analyses.
Result Summary	Opens the Summaries view.
Result Filters	Opens the Result Filters view for reducing the number of rows displayed in selected result tables.
Chromatograms	Opens the Chromatograms view for viewing chromatogram plots.

Menu command	Description	
Mass Spectrum	Opens the Mass Spectrum view for viewing a spectral tree and th spectrum scans.	
Trend Chart	Opens the Trend Chart view for setting up a box-and-whisker (Box Whisker selection) chart or a trendline chart.	
Isotopologues Distribution Chart	Opens the Isotopologues Chart view. Available only when the active result file includes results from the Analyze Labeled Compounds node.	
Show Mass Defect Plot	Opens the Mass Defect Plot view for examining the relationship between the mass defects of the detected or expected compounds and their molecular weights.	
Result Charts	Opens the Result Charts view for setting up data graphs, such as scatter plots, histogram charts, bar charts, and pie charts. Use these views to visualize the data.	
Descriptive Statistics	Opens the Descriptive Statistics view for viewing a box-and-whisker plot of all the compounds in the Compounds or Expected Compounds tables for the selected sample groups.	
Differential Analysis	Opens the Differential Analysis view for viewing volcano plots and running differential analyses.	
Principal Component Analysis	Opens the Principal Component Analysis view for evaluating multivariate data.	
Partial Least Squares Discriminant Analysis	Opens the Partial Least Squares Discriminant Analysis view.	
Hierarchical Cluster Analysis	Opens the Hierarchical Cluster Analysis view.	
Retention Time Corrections	Opens the Corrected Retention Times view. Available only when the active result file includes data from multiple input files. To view the retention time correction cu for one or more input files, select the input files of interest in Input Files table.	
Compound Area Corrections	Opens the Compound Area Corrections view. Available only when the active result file includes data from Quality Control samples.	
Metabolika Pathways	Opens the Metabolika Pathways view for viewing the Metabolika pathways that are mapped to the compounds data.	

Table 4. Application menu bar (Sheet 4 of 6)

Menu command	Description	
BioCyc Pathways	Opens the BioCyc Pathways view for viewing the BioCyc pathways that are mapped to the compounds data.	
	Available only when the active result file includes mapped BioCyc pathways.	
KEGG Pathways	Opens the KEGG Pathways view for viewing the KEGG [™] pathways that are mapped to the compounds data.	
	Available only when the active result file includes mapped KEGG pathways.	
mzLogic Analysis	Opens the mzLogic Analysis view.	
FISh Scoring Queue	Opens the FISh Scoring Queue view.	
Window menu		
Use these command to a	apply, save, manage, or reset layouts.	
Apply Layout	Displays the layouts list for selecting a layout.	
Save Layout	Opens the Save Layout dialog box for naming the current layout.	
Manage Layouts	Opens the Manage Result File Layout dialog box for renaming or deleting layouts.	
Reset Layout	Closes and reopens the active result file to reset its layout.	
Help menu		
These commands are alw	ways available.	
Compound Discoverer Help	Opens the Compound Discoverer Help, which is a compiled Help file with Contents, Index, and Search tabs. The Help contains context-sensitive topics; that is, pressing F1 opens the Help topic that corresponds to the current area of the application.	
How to Use the Help	Opens the Compound Discoverer Help to the Using This Help topic.	
Glossary	Opens the Compound Discoverer Help to the table of contents for the glossary.	
Compound Discoverer Support Manuals	Displays links to the Compound Discoverer User Guide, Compound Discoverer Reporting Quick Start, Compound Discoverer tutorials, and Release Notes.	
License Manager	Opens the License Manager page where you can activate your Compound Discoverer license or scan for missing application features.	

Table 4. Application menu bar (Sheet 5 of 6)

Menu command	Description
Communication Tests	Opens the Communication Tests dialog box where you can run tests to check whether your data processing computer can connect to the external BioCyc Pathways, KEGG Pathways, mzCloud, and ChemSpider databases.
Create Bug Report	Creates a report of your computer's configuration and stores it as a Compound Discoverer Bug Report (<i>timestamp</i>).zip on your computer desktop.
	To report software errors to Thermo Fisher Scientific, send a detailed error description with screen shots and attach this bug report to the email.
Configuration	Opens the Configuration page for setting global options, such as the maximum number of parallel analyses, the study management setting for automatically saving studies, the color maps, the mzCloud search settings, and your user credentials for the BioCyc mapping feature.
About	Opens the Compound Discoverer dialog box for viewing lists of the installed components and processing workflow nodes.

Table 4. Application menu bar (Sheet 6 of 6)

Application Toolbar

Figure 4 shows the application toolbar.

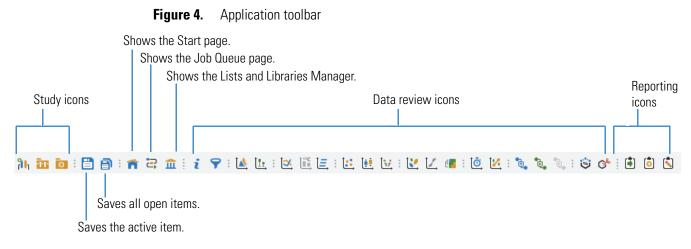


Table 5 describes the icons in the application toolbar from left to right.

Table 5.	Toolbar icons (Sheet 1 of 3)
----------	------------------------------

lcon	Description		
Study ico	ns		
ř.	Opens the New Study and Analysis Wizard that takes you through the proc of specifying the studies folder for all your study subfolders, creating a new study, and starting a new analysis.		
æ	Opens the Open Study dialog box for selecting the current version or a previous version of an existing study.		
0	Opens the Open Result File dialog box for selecting the current version or a previous version of a result file.		
General i	cons		
	Saves the currently active item, such as a study or result file.		
ð	Saves all the open pages, such as the study pages and the result pages.		
n	Opens the Start Page if it is not already open and makes it the active page.		
2	Opens the Job Queue page if it is not already open and makes it the active page.		
血	Opens the Lists and Libraries Manager where you can select to open one of th following lists or libraries:		
	• Expected Compounds—For viewing or modifying a list of compounds		
	• Transformations—For viewing or modifying a list of transformations		
	• Ion Definitions—For viewing or modifying a list of ion definitions		
	• Adducts—For viewing or modifying a list of adducts		
	• Mass Lists—For creating, editing, importing, exporting, or deleting mass lists		
	• Spectral Libraries—For creating, editing, importing, exporting, or deletin mzVault database files		
	 Metabolika Pathways—For adding, editing, importing, exporting, or deleting Metabolika pathways. 		
	 Compound Classes—For creating, editing, importing, exporting, or deleting compound class libraries of fragment structures 		

Table 5.Toolbar icons (Sheet 2 of 3)

lcon	Description		
Data review	/ icons		

Available when a result file is the current page in the application window. When the respective view is open, brings the view to the forefront or makes the view the active view.

1	
i	Opens the Summaries view as a docked view. Available when the Summaries view is closed.
9	Opens the Result Filters view as a floating window.
?	Opens the Chromatograms view as a docked view.
ţ ļ	Opens the Mass Spectrum view as a docked view.
î 🏹	Opens the Trend Chart view as a docked view.
Ĵ∎ [™] c	Opens the Isotopologues Distribution Chart view as a docked view.
••	Available when the opened result file includes an analysis of labeled compounds.
î <i></i> ,	Opens the Mass Defect Plot view as a docked view.
Î.	Opens the Results Chart view as a floating window.
<u> </u>	The Results Chart view includes the following pages: Scatter Chart, Histogram Chart, Bar Chart, and Pie Chart.
Ĵ∎ !	Opens the Descriptive Statistics view as a docked view.
ÎV,	Opens the Differential Analysis view as a docked view.
Ì¥,	Opens the Principal Component Analysis view as a docked view.
	Opens the Partial Least Squares–Discriminant Analysis (PLS–DA) view as a docked view.
(Opens the Hierarchical Cluster Analysis view as a docked view.
<u>(م)</u>	Opens the Retention Time Corrections view as a docked view. Available when the result file includes more than one input file.

lcon	Description	
Î::/;	Opens the Compounds Area Correction view as a docked view.	
1	Opens the Metabolika Pathways view as a docked view. Available when the result file includes mapped BioCyc pathways.	
Ð	Opens the BioCyc Pathways view as a docked view. Available when the result file includes mapped BioCyc pathways.	
B	Opens the KEGG Pathways view as a docked view. Available when the result file includes compounds mapped to KEGG pathways.	
	Selecting an item in the KEGG Pathways result table opens the reference pathway for the item in the KEGG Pathway view.	
ð	Opens the mzLogic Analysis view as a docked view.	
ß	Opens the FISh Annotations Queue view as a docked view.	
Reporting	icone	

Table 5. Toolbar icons (Sheet 3 o	f 3)
-----------------------------------	------

Available when a result file (not necessarily the current page) is open in the application window.

.	Opens the Open Report Design Template dialog box for selecting a report template (.cdReportTemplate).
ी	Opens the Customize Report dialog box for setting up a custom report template.
(Opens the Open Report Design Template dialog box for selecting a report template (.cdReportTemplate).

Understanding Processing Workflows

The application uses a node-based method to create processing workflows for the analysis of raw data files. You can create your own custom processing workflows (cdProcessingWF), use one of the many defined processing workflow templates provided with the application, or customize one of the existing processing workflow templates.

Note The parameter settings for the workflow nodes in the defined processing workflow templates are not necessarily the same as the default settings for the workflow nodes themselves, as Thermo Fisher Scientific has optimized the parameter settings in the defined processing workflow templates by the area of study (vertical market).

These topics describe the defined processing workflow template and how the application processes data to find expected compounds for a targeted analysis or detect and identify unknown compounds for an untargeted analysis:

- Targeted Workflows for Expected Compounds
- Untargeted Workflows for Identifying Unknown Compounds
- Marking Background Compounds in an Untargeted Analysis
- Nomenclature for the Defined Processing Workflows
- The Defined Processing Workflows

Related Topics

- Workflow Nodes
- Creating and Editing Processing Workflows

Targeted Workflows for Expected Compounds

Use the Expected Compounds workflow nodes to run an analysis that targets known analytes—for example, the metabolites of a specific drug.

Figure 5 shows the workflow tree in the following processing workflow file:

MetID w Stats Expected w FISh Scoring and Background Removal.cdProcessingWF

This processing workflow uses the Generate Expected Compounds and Find Expected Compounds nodes to find expected parent¹ compounds and their dealkylation, dearylation, and transformation products. FISh Scoring is applied to explain the fragments in the fragmentation spectra based on in silico fragment prediction of the parent and dealkylation/dearylation products.

¹ A parent compound is the initial compound in a reaction or metabolic pathway.

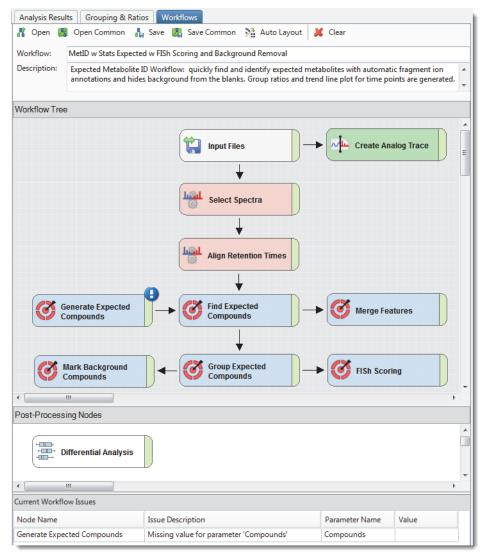


Figure 5. Processing workflow for a metabolism study

During this targeted analysis, the following processes occur:

- 1. The Input Files Node sends the file names and location of the input files to the connected nodes (typically the Select Spectra node, Create Analog Trace node, or both of these nodes).
- 2. The Select Spectra Node filters the MS scan data.

IMPORTANT For a targeted analysis, which uses the Find Expected Compounds node, do not filter out all the full (MS1) scans, as this node requires the full scan data. When the processing workflow includes the FISh Scoring node, do not filter out the fragmentation scans, as this node uses them to provide a confirmation score for the expected precursor ions.

3. The Align Retention Times Node chromatographically aligns features across the input files in a sample set by using the specified alignment algorithm.

IMPORTANT The alignment algorithm looks for matching features (chromatographic peaks with the same $m/z \times RT$ dimensions) in the input files. The alignment algorithm can align input files that include polarity switching data; it cannot align input files that include only positive polarity scans with input files that include only negative polarity scans.

- 4. The Generate Expected Compounds Node creates a list of expected compound ions by using the following user-specified parameter settings:
 - Parent compound or compounds
 - Number of dealkylation and dearylation steps
 - Number and type of transformation steps
 - List of possible adduct ions

The Generate Expected Compounds node passes the following information to the Find Expected Compounds node for each expected compound:

- Parent compound or compounds
- Elemental composition of the expected neutral compound
- Molecular weight of the expected neutral compound
- Whether the expected compound is a product of a dealkylation step
- Transformations required to produce the expected compound
- Composition change between the parent and expected compound

It also passes the following information for each expected compound ion:

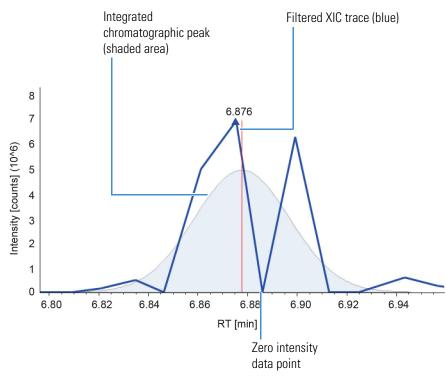
- Charge
- Theoretical *m/z* value

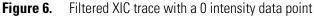
Note The Generate Expected Compounds node does not send information to a result table. It sends information to the Find Expected Compounds node, which requires input from at least one Generate Expected Compounds node. The Filter By Mass Defect node can also process the input from one or more Generate Expected Compounds nodes.

- 5. The Find Expected Compounds Node performs the following steps by using the full (MS1) scan data from the Align Retention Times node and the information from the Generate Expected Compounds node (or multiple Generate Expected Compounds nodes):
 - a. Creates a set of mass tolerance and intensity tolerance fit parameters for each expected compound ion (m/z value) by using the theoretical m/z value for the ion, the user-specified mass tolerance, the theoretical isotope pattern for the ion, and the user-specified intensity tolerance for the isotopic ions.
 - b. For each expected compound ion, the Find Expected Compounds node does the following:
 - i. Checks each full (MS1) scan that passes through the connected data processing node for centroids that match the mass and intensity tolerance rectangles. The pattern search (set of rectangles) looks for the base peak (most intense centroid) of the pattern first.

In most cases, the base peak is the A0 centroid for the monoisotopic ion. But in some cases, for example, in compounds that contain two bromine atoms or four chlorine atoms, the isotopic peaks have a higher intensity than the monoisotopic peak.

- ii. Draws a filtered XIC trace by summing the centroids found for each data point. If a data point does not contain the user-specified number of matching isotopes and the theoretical intensity of the missing isotope was above the noise threshold, the node assigns a 0 intensity value to the data point (Figure 6).
- c. For each expected compound found, the Find Expected Compounds node creates a summed trace by summing the XIC traces of the associated expected compound ions.
- d. Detects and integrates the chromatographic peaks in each XIC trace. Does not report chromatographic peaks with an apex peak height that is below the user-specified minimum (chromatographic) peak intensity. If the average peak width of the peaks in the processed retention time range is greater than the setting for the Average Peak Width parameter, the Find Expected Compounds node rejects all of the chromatographic peaks.





Note A chromatographic peak with a 0 intensity value indicates one of the following:

• The user-specified mass tolerance, intensity tolerance, or number of required isotopes for the Find Expected Compounds node is not suitable for this sample set. Modify these parameter settings and rerun the analysis.

-or-

• The chromatographic peak does not represent the presence of an expected compound. Ignore this chromatographic peak.

The Find Expected Compounds node passes the following information to the Expected Compounds per File table for each expected compound that it finds in an input file:

• RT (min)—Retention time of the chromatographic peak apex.

Note If the node finds more than one adduct ion for an expected compound, the chromatogram is a summed trace.

- Best SFit [%]—Best spectral fit score between the measured and expected isotope patterns for the expected compound ions. When the node finds only one adduct ion, the best spectral score is equal to the score for the adduct ion that it found.
- Max # MI—Maximum number of matched isotopes for any of the adduct ions. When the node finds only one adduct ion, the maximum number of matched isotopes is equal to that of the adduct ion that it found.

- #Adducts—Number of adduct ions that it found.
- Area—Total area of the chromatographic peaks for the found adduct ions of the expected compound. Peak areas are reported in counts × seconds.
- Parent Area%—Relative area of the chromatographic peak for the expected compound as compared to the total area of all found peaks for the expected compound.

Note The parent area is the chromatographic peak area of the expected compound, rather than the area of the compound listed in the Parent Compound column. The compounds listed in the Parent Compound column are the library compounds that you selected in the Generate Expected Compounds node.

6. The Group Expected Compounds Node groups the chromatographic peaks by their molecular weight × retention time (MW × RT) dimensions across the input file set and creates the Expected Compounds table.²

This node also selects the best hit ions for each compound across the input file set:

- Selects the best ion and related MS1 scan for each compound as the one with the highest resolution and the highest intensity for the preferred ion. When you open a result file, the mass spectrum view displays the MS1 scan for the best ion across the input file set.
- Selects the best fragmentation data by using the user-specified preferred precursor ion with the highest intensity that has data-dependent MS2 scans.

Note You use the Preferred Ions parameter in the Group Expected Compounds node to specify the preferred adduct ions.

- 7. When the input file set includes Blank samples (Sample Type: Blank), the Mark Background Compounds node compares the peak areas of the compounds (same parent compound, molecular weight, and retention time) that are found in both the Blank samples and the non-blank samples, and labels these compounds as Background Compounds if their peak areas do not meet the specified threshold.
- 8. When the spectrum data in the input files includes fragmentation scans (MS2, MS3, MSn), the FISh Scoring node compares the best fragmentation scan for an expected compound across the input file set to the expected structures. The scoring process can add considerable processing time.
- 9. When the processing workflow includes both the Find Expected Compounds node and the Detect Compounds node, the Merge Features node consolidates the chromatographic peaks that these nodes find in the main Merged Features table. The consolidation is based on the $m/z \times RT$ dimensions of the features.

² For result files generated in Compound Discoverer 2.0, the Group Expected Compounds node also added the Best Compounds column to the Expected Compounds per File table. The Workflow page of the Result Summary view displays the Best Compound criteria for updated result files.

Note The Merge Features node also creates the Merged Features table that is related to the Manual Peaks table. When you manually integrate a chromatographic peak for a specialized trace such as a UV trace, you can compare the selected peak in the Manual Peaks table to the peaks in the related Merged Features table. The application populates the related Merged Features table with the chromatographic peaks that fall within the retention time window specified in the Merge Features node—that is, the application populates the Merged Features table that is related to the Manual Peaks table with chromatographic peaks that have a similar retention time to the selected manual peak.

10. When the analysis includes group ratios, the Differential Analysis node runs the differential analysis.

After the analysis finishes, you can open the result file.

Figure 7 shows a schematic of the main and related result tables for the targeted workflow shown in Figure 5. The Structure Proposals table is empty until you populate it. For information about adding structure proposals or editing the compound annotations, see "Editing Compound Annotations" on page 259.

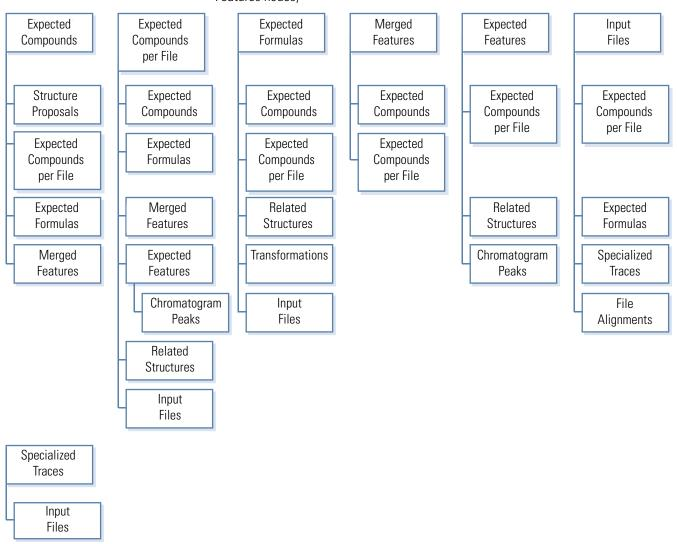


Figure 7. Result tables for a basic targeted workflow (with the Create Analog Traces and Merge Features nodes)

Untargeted Workflows for Identifying Unknown Compounds

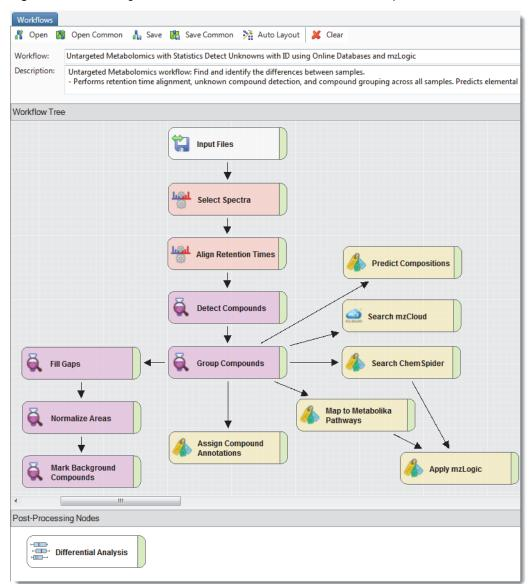
Figure 8 shows the workflow tree for the metabolomics processing workflow that is provided with the application:

Untargeted. Metabolomics with Statistics Detect Unknowns with ID using Online Databases and mzLogic.cdProcessingWF

This processing workflow uses the Detect Compounds node to find chromatographic peaks for unknown compounds (MW \times RT) and the Predict Compositions node to determine the possible elemental compositions of the unknown compounds. It also determines the possible identity of the unknown compounds as follows:

- The Search ChemSpider node searches selected databases of MS1 scans by using the MW or predicted formulas when available.
- The Search mzCloud node searches the mzCloud database of fragmentation scans by using the MW or predicted formulas when available.
- The Map to Metabolika Pathways searches the pathways for the detected compounds. When the application finds a matching compound, it maps the pathway for ease of viewing.

The Assign Compound Annotations node assigns the following annotations to the detected compounds: Name, Formula, and Structure. The Apply mzLogic node combines mzCloud similarity searching (MS2 and MSn) with structure similarity matching to rank putative database results.





During this untargeted analysis, the following processes occur:

- 1. The Input Files node sends the file names and location of the input files to the Select Spectra node.
- 2. The Select Spectra node filters the MS scan data.
- 3. The Align Retention Times node chromatographically aligns the input files in a sample set by using the specified alignment algorithm.
- 4. The Detect Compounds node detects XIC traces in the full (MS1) scans by using the parameter settings for the mass tolerance and intensity threshold, groups isotopes and adducts by using the user-specified ions and base ions lists, and reports the unknown compounds (MW×RT) by occurrence in the Compounds per File table.

IMPORTANT Make sure that the Preferred Ions list for the Group Compounds node includes the selections in the Base Ions list for the Detect Compounds node.

5. The Group Compounds node uses the specified mass and RT tolerances to group chromatographic peaks with the same MW×RT values in the Compounds table. It then sends the best fragmentation data across the input files to the Search mzCloud node and Predict Compositions node.

This node also selects the best hit ions for each compound across the input file set:

- Selects the best ion and related MS1 scan for each compound as the one with the highest resolution and the highest intensity for the preferred ion. When you open a result file, the mass spectrum view displays the MS1 scan for the best ion across the input file set.
- Selects the best fragmentation data by using the user-specified preferred precursor ion with the highest intensity that has data-dependent MS2 scans.

The Predict Compositions node and the search nodes use the best hit ions.

Note You use the Preferred Ions parameter in the Group Compounds node to specify the preferred adducts.

- 6. The Search mzCloud node searches the mzCloud database for matching and similar fragmentation spectra.
- 7. The Predict Compositions node predicts the elemental compositions of the unknown compounds.
- 8. The Map to Metabolika Pathways node searches the Metabolika pathways for matching compounds.
- 9. The Search ChemSpider node searches the ChemSpider database for matching compounds.
- 10. The Fill Gaps node fills in missing peaks or peaks below the detection threshold (specified in the Detect Compounds node) for subsequent statistical analysis.
- 11. The Mark Background Compounds node determines the background compounds in the Blank samples and labels these compounds as background compounds.
- 12. The Differential Analysis node runs a differential analysis on the defined sample ratios and calculates the p-values.
- 13. The Assign Compound Annotations node assigns and compares the annotations provided by the Predict Compositions, Search ChemSpider, Search mzCloud, and Search Mass Lists nodes.

14. The Apply mzLogic node runs a forward search and a reverse search using the mzCloud service. For compounds that have available MS2 scans, it scores all the structure candidates (or the specified maximum number of candidates) from the attached input nodes. Adds the following columns to the result tables: #Similarity Results (Compounds table) and mzLogic Score (search result tables).

For information about the result tables that this processing workflow generates, see "Compound Detection Result Tables" on page 401.

Figure 9 shows a schematic of the main and related tables for a basic untargeted workflow.

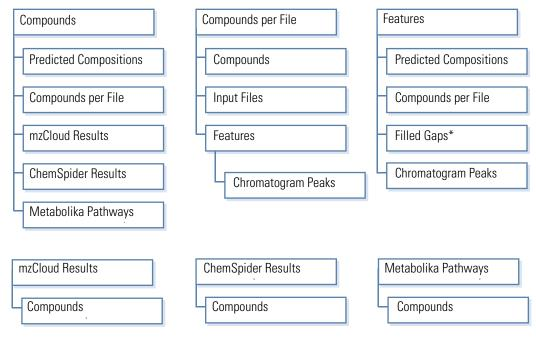


Figure 9. Result tables for an untargeted workflow

*By default, the Filled Gaps table is hidden.

Marking Background Compounds in an Untargeted Analysis

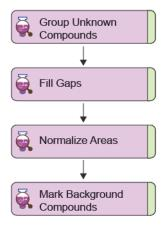
This topic describes how to make the appropriate node connections for an untargeted processing workflow that includes two or all three of these workflow nodes: Fill Gaps, Mark Background Compounds, and Normalize Areas. All three of these nodes process data across the input file set.

The Fill Gaps and Normalize Areas nodes automatically make the appropriate connections:

- The Fill Gaps node only takes input from the Group Compounds node. When you add the Fill Gaps node to a processing workflow that has a Group Compounds node, the Group Compounds node automatically connects to it.
- The Normalize Areas node only takes input from the Fill Gaps node. When you add the Normalize Areas node to a processing workflow that has a Fill Gaps node, the Fill Gaps node automatically connects to it.

The Mark Background Compounds node can take input from any of these three nodes— Group Compounds, Fill Gaps, or Normalize Areas. However, always add the Mark Background Compounds node to the end of the processing workflow.

When an untargeted processing workflow includes the Fill Gaps, Normalize Areas, and Mark Background Compounds nodes, the correct connection order is as follows: Group Compounds > Fill Gaps > Normalize Areas > Mark Background Compounds.



Related Topics

- Fill Gaps Node
- Normalize Areas Node
- Mark Background Compounds Node–Unknown Compounds

Nomenclature for the Defined Processing Workflows

When you open a defined processing workflow on the Workflows page of an analysis, descriptive text appears in the Description box. Read and understand the description before you use the workflow. The Description box does not enlarge to fit the text, so you must scroll down to read the complete description. Or, you can click the Description box, press CTRL + A to copy the text to the Clipboard, and paste the text into Notepad.

The file names for the processing workflows include the following descriptive text:

• Expected—The workflow runs a targeted analysis with the Generate Expected Compounds and Find Expected Compounds nodes.

IMPORTANT You must customize the targeted workflows by selecting the targeted compounds in the Generate Expected Compounds node. Before you can select the compounds for the node, you must add the compounds to the Expected Compounds library.

In the defined processing workflows for targeted analyses, the Generate Expected Compounds node is set up to generate a mass list for the following adduct ions: [M+H]+1 and [M–H]–1. For best results, make the appropriate selections for your analysis from the Ions list in the Generate Expected Compounds node.

• Detect Unknowns, Unknown, or Untargeted—The workflow runs an untargeted analysis with the Detect Compounds node.

Tip Mobile phase additives can have a significant effect on the base ions (adduct ions with the highest intensity) in the full scan data for an LC/MS experiment. To avoid misinterpreting the isotopic ion clusters, make sure that the Base Ions list in the Detect Compounds node includes the predominant adduct ions. For example, if the mobile phase contains a significant amount of ammonium acetate, consider adding the ammonium adduct, [M+NH4]+1, to the list.

The processing workflows in the E and L folder include [M+NH4]+1 in the Base Ions list.

• Online Databases—The workflow searches the mzCloud and ChemSpider databases.

Tip In the Search ChemSpider node, select the appropriate databases.

• Local Database—The workflow searches your local mzVault and mass list files. For metabolomics templates, the local databases include the local Metabolika pathways files.

Tip In the Search mzVault and Search Mass Lists nodes, select the appropriate files. The application includes four mzVault library files and four mass lists.

- mzVault libraries:
 - mzVault Reference May 2019 (snapshot of the mzCloud spectral database, MS2 data only)
 - mzVault Autoprocessed May 2019 (snapshot of the mzCloud spectral database, MS2 data only)
 - Bamba Lab 34 Lipid Mediators Library Stepped NCE 10 30 45
 - Bamba Lab 598 Polar Metabolites Stepped NCE 10 30 45
- mass lists
 - Arita Lab 6549 Flavonoid Structure Database
 - EFS HRAM Compound Database
 - Endogenous Metabolites Database 4400 Compounds
 - Extractables and Leachables HRAM Compound Database
- FISh Scoring—The workflow includes the FISh Scoring Node, a structural confidence scoring and annotation tool for comparing the predicted fragments of expected compounds to the experimental fragmentation scans. This node adds a significant amount of processing time.
- Stats or Statistics—The workflow includes the Differential Analysis node. If you submit an analysis that does not include ratios, a warning message appears. If you do not want to run a differential analysis, you can ignore the warning and submit the run.
- Compound Class Scoring—The workflow includes the Compound Class Scoring Node.

IMPORTANT Before you can select the compounds classes for the node, you must add the fragment lists to the Compound Classes library.

- MMDF—The workflow includes the Filter By Mass Defect Node.
- Background Removal—The targeted workflow includes the Mark Background Compounds node for filtering out expected compounds that are also found in the blank samples.

Note All the untargeted workflows include the Mark Background Compounds node for filtering compounds detected in the blank samples.

• Pattern Scoring—The workflow includes the Pattern Scoring node.

The Defined Processing Workflows

Tip To access these workflows from the Workflows page of an analysis, click **Open Common** from the command bar, open the Workflow Templates folder, open the folder for the applicable vertical market, and then select a template.

The processing workflows installed with the application are as follows:

- DegradantID folder
 - Degradants Related and Unknown ID w Database Searches
 - Degradants Unknown ID w Pattern Trace and Pattern Scoring
 - Degradants w Stats Related and Unknown ID w Database Searches
- EandL folder
 - E and L Expected w FISh Scoring
 - E and L Unknown ID with Online and Local Database Searches
 - E and L w Stats Unknown ID w Online and Local Database Searches
- Environmental folder
 - Environmental Expected w Transformation and FISh Scoring
 - Environmental Unknown ID w Online and Local Database Searches
 - Environmental w Stats Unknown ID w Online and Local Database Searches
- Food Research folder
 - Food Research Expected w FISh Scoring
 - Food Research Unknown ID w Online and Local Database Searches
 - Food Research w Stats Unknown ID w Online and Local Database Searches
- Forensics folder
 - Forensics Expected w FISh Scoring
 - Forensics Unknown ID w Compound Class Scoring and Database Searches
 - Forensics w Stats Unknown ID w Compound Class Scoring and Database Searches
- ImpurityID folder
 - Impurity ID Related and Unknown
 - Impurity ID w Stats Related and Unknown

- Metabolomics folder
 - Max ID Detect Unknowns with ID Using Online Database Searches Single Sample
 - Untargeted Metabolomics Quick Detection Unknowns No ID
 - Untargeted Metabolomics using Online Databases, mzLogic, and Molecular Networks
 - Untargeted Metabolomics with Statistics Detect Unknowns with ID using Local Databases
 - Untargeted Metabolomics with Statistics Detect Unknowns with ID using Online Databases
 - Untargeted Metabolomics with Statistics Detect Unknowns with ID using Online Databases and mzLogic
- MetID folder
 - MetID Generate Inclusion List For Acquisition Pos Mode
 - MetID Generate Inclusion List For Acquisition Neg Mode
 - MetID Pattern Scoring with Background Removal
 - MetID w Stats Expected and Unknown w Background Removal
 - MetID w Stats Expected and Unknown w MMDF and Background Removal
 - MetID w Stats Expected w Background Removal
 - MetID w Stats Expected w FISh Scoring and Background Removal
- NaturalProduct
 - Natural Product Unknown ID w Online and Local Database Searches
 - Natural Product Unknown ID w Stats Online and Local Database Searches
- PolymerID
 - Unknown Polymer ID w Stats Online and Local Database Searches
 - Unknown Polymer ID with Online and Local Database Searches
- Stable Isotope Labeling folder
 - Stable Isotope Labeling w Metabolika Pathways and ID using Local Databases
 - Stable Isotope Labeling w Metabolika Pathways and ID using Online Databases

Best Scans for Composition Prediction and Spectral Matching

For each set of raw data files that you submit for processing, the application can select the best MS1 scan and the best MS2 scan for each compound from different input files. This selection is done within the compound consolidation nodes—that is, by the Group Compounds node and the Group Expected Compounds node.

For information about how the application determines and uses the best scans, see these topics:

- Best MS1 Scan for Isotope Pattern Matching
- Best MS2 Scan for Fragments Matching and Spectral Comparison

Best MS1 Scan for Isotope Pattern Matching

To determine the best MS1 scan to use for a compound across a set of input files, the consolidation nodes select the MS1 scan with the highest resolution and retention time closest to the compound's peak apex.

The application sends the isotope pattern information from the best MS1 scan to the following workflow nodes for further processing:

- To the Predict Compositions node for isotope pattern matching
- To the Apply Spectral Distance node for calculating the spectral distance between the assigned elemental composition and the experimental isotope pattern

Best MS2 Scan for Fragments Matching and Spectral Comparison

To identify compounds by fragmentation data or to determine the spectral similarity between the fragmentation scans for two different compounds, the application uses the "best MS2 scan" for each compound.

To determine which MSn scans to attach to a compound, the application does the following:

- 1. Searches for MS2 scans for the precursor ion within the RT range of the compound's peak apex ± FWHM.
- 2. If it finds no MS2 scans for the precursor ion within this range, it searches for scans within the start and end points of the compound's chromatographic peak, as determined by the Parameterless Peak Detection (PPD) algorithm.
- 3. To determine the best MS2 scan, it does the following:
 - a. Selects the MS2 scan for the preferred ion.
 - b. If it finds multiple spectrum trees for the preferred ion, it uses the MS2 scan from the MS1 scan with the highest precursor intensity.

For each compound, the Group Expected Compounds and the Group Compounds nodes send all the MS2 scans associated with that compound to the following nodes:

- Search mzCloud node
- Search mzVault node

For each compound, the Group Compounds node sends the best MS2 scans to the Apply mzLogic node.

Quality Control Samples for Batch Normalization

When acquiring raw data files for a large sample set, use pooled quality control samples to compensate for time-dependent batch effects. For information about batch effects, refer to the following article: *Procedures for large-scale metabolic profiling of serum and plasma using gas chromatography and liquid chromatography coupled to mass spectrometry*.³

To create a pooled quality control sample, combine a small aliquot from each sample in the processing batch (input file set to be processed to create one result file). When setting up an acquisition sequence, bracket the unknown samples by injecting this pooled quality control sample at regular intervals.

When the processing workflow includes the Normalize Areas Node and the input file set includes QC samples, the application uses the QC samples to create a regression curve of area versus acquisition time for each detected compound.

The application does not create a regression curve for a particular compound and does not correct the areas in the non-QC samples unless all three of these conditions are met:

- It detects the compound in the user-specified minimum percentage of the QC samples.
- The relative standard deviation of the detected peak areas for the compound in the QC samples does not exceed the user-specified threshold.

Note The application checks the RSD of the peak areas both before and after applying a correction for the relative acquisition time. If the RSD for the corrected peak areas differs by more than 75% of the user-specified threshold, the application does not create a corrected-area regression curve for the compound.

• The number of samples acquired between the QC samples does not exceed the user-specified number.

³ Dunn, W.B.; Broadhurst, D.; Begley, P.; Zelena, E.; Francis-McIntyre, S.; Anderson, N.; Brown, M.; Knowles, J.D.; Halsall, A.; Haselden, J.N.; Nicholls, A.W.; Wilson, I.D.; Kell, D.B.; Goodacre, R. Human Serum Metabolome (HUSERMET) Consortium. *Nat Protoc.* **2011**, 6(7), 1060-83.

You can view the results of the batch normalization process in the Compounds result table and the Compound Area Corrections view. The Compounds table displays the corrected compound areas. The Compound Area Corrections view shows the effect of the signal correction on the absolute compound areas. If the application excluded the selected compound during the QC signal correction step, the Compound Area Corrections view is empty.

Stable Isotope Labeling

Stable isotope labeling experiments are fundamental to understanding metabolic pathways and the turnover rate (flux) for the molecules in these pathways. Where traditional approaches are usually limited to targeted analyses, which detect and quantify known compounds and their labeled downstream metabolites, the Compound Discoverer application makes full use of high-resolution Orbitrap[™] data and offers state-of-the-art analyses for unknown labeled compounds.

Using an unlabeled reference sample, the application detects unknown compounds above a specified minimum intensity threshold, determines their elemental composition and identity, and then determines the labeled counterparts (isotopologues) of these compounds in the samples marked as labeled.

You can use any isotopic label in your experiments; however, when you use a label other than carbon-13, you must specify the labeled element in the processing workflow.

The application reports the isotopologues and the fractional label incorporation (exchange rate) for each compound. You can overlay the exchange rate as well as other statistical data onto pathways using Metabolika[™], which is directly integrated into the Compound Discoverer application.

To acquire and process stable isotope labeled data, do the following:

- 1. Acquire LC/MS/MS data for a set of samples where at least one sample corresponds to the unlabeled state of the system.
- 2. Select the following sample types:
 - For unlabeled samples, select **Sample**, **Identification Only**, **Quality Control**, or **Blank** as appropriate. You must select Sample (or Control or Standard) for at least one sample.

Note Currently, the application treats Sample, Control, and Standard samples the same way. This functionality is subject to change in future releases.

- For labeled samples, select Labeled.
- 3. Select one of the processing workflows in the following folder:

Common Templates > Workflow Templates > Stable Isotope Labeling

4. When using a label other than carbon-13, customize the parameter settings for the Analyze Labeled Compounds node.

During data processing, the application does the following:

- 1. Detects and identifies the compounds in the unlabeled samples.
- 2. Submits the compounds with assigned formulas to the Analyze Labeled Compounds node which does the following:
 - a. Generates isotopologues for each detected compound by replacing one occurrence of an atom at a time by its specified isotope. The compound's formula or the user-specified limit defines the number of exchangeable atoms, whichever is smaller.
 - b. Simulates the isotope pattern for each isotopologue by using its formula and the spectral resolution of the raw data.
 - c. Consolidates the isotopologue patterns for each compound to get the final set of expected masses, considering the specified mass tolerance and the spectral resolution of the raw data.
 - d. Generates an XIC trace for each expected mass and detects the chromatographic peaks.
 - e. Deconvolves the chromatographic peaks to determine the relative amount of each isotopologue.
 - f. Flags compounds containing contaminating masses in unlabeled samples, as well as unusual isotopologue distributions and insufficient pattern fits.

FISh Scoring for Targeted Compounds and Proposed Structures

Note In the current version of the application, the FISh scoring algorithm uses all the fragmentation scans (in the spectrum tree for a compound), compared to earlier versions of the application that only used MS2 scans. Therefore, if you reprocess data sets that you already processed with an earlier version of the application, and these data sets include MSn scans (where n > 2), the FISh coverage scores might be lower.

For targeted compounds, you can add FISh scoring to the automated analysis, and for unknown compounds, you can apply FISh scoring to your proposed structures.

The FISh scoring algorithm attempts to match the fragment structures in a list of expected fragments to the centroids in the fragmentation scans of the precursor ions.

When a precursor ion scan is followed by only one fragmentation scan, the node calculates the FISh coverage score as follows:

FISh coverage score = $\frac{\# \text{ matched centroids}}{\# \text{ used centroids}} \times 100$

where:

matched centroids represents the number of matched centroids.

used (matched + unmatched) centroids represents the number of centroids in the fragmentation scan that are above the user-specified signal-to-noise threshold. The algorithm skips centroids below the user-specified signal-to-noise threshold.

When a precursor scan is followed by more than one fragmentation scan, the node calculates a composite score as follows:

FISh coverage score =
$$\frac{(\sum_{\text{per all scans}} \# \text{ matched centroids})}{(\sum_{\text{per all scans}} \# \text{ used centroids})} \times 100$$

The FISh scoring algorithm annotates the centroids in the fragmentation scans with the matching fragment structures. It also provides a FISh Coverage score for data-dependent scans in the Mass Spectrum view legend and a FISh Coverage score in the Expected Compounds table.

Calculating the Mass Defect of an Elemental Composition

In the Compound Discoverer application, you can use the mass defect of an elemental composition as follows:

- Filter the spectral data during processing to keep or remove expected and detected compounds by their mass defect.
- Calculate the mass defect of each detected compound by using multiple calculation methods during processing, and then sort the resulting table of detected compounds by their mass defect.
- View a plot of the mass defects for the detected compounds versus their molecular weights to visualize the similarity between the compounds.

The application uses the following formulas to calculate the mass defect of an elemental composition.

Mass defect type	Formula	
Fractional mass	exact mass – floor of the exact mass	
	where:	
	exact mass = monoisotopic mass of the elemental composition	
Standard mass defect	exact mass – nominal mass	
	where:	
	nominal mass = integer mass of the elemental composition	
	The application calculates the integer mass by using the selected rounding function:	
	• Floor rounds down	
	• Ceiling rounds up	
	• Round rounds to the nearest integer value	
Relative mass defect	1e ⁶ × (exact mass – nominal mass)/exact mass	
Kendrick mass defect	Kendrick mass – nominal Kendrick mass	
	where:	
	Kendrick mass = $a \times (b/c)$	
	a = exact mass of the elemental composition	
	b = nominal mass of the Kendrick formula	
	c = exact mass of the Kendrick formula	

Table 6. Mass defect types

Related Topics

- Filter By Mass Defect Node
- Calculate Mass Defect Node
- Working with the Mass Defect Plot View

Using mzLogic to Score Candidates for Unknown Compounds

There are two ways to run the mzLogic scoring algorithm:

- Running an Untargeted Analysis that Includes the mzLogic Node
- Running an mzLogic Analysis from the mzLogic Analysis View

Running an Untargeted Analysis that Includes the mzLogic Node

Processing workflow templates that include the following terms—Database Searches or mzLogic—in their file names include the mzLogic node and one or more of the compound identification nodes and pathway mapping nodes.

When an mzCloud search yields no identity matches for an unknown compound, the mzLogic algorithm provides a ranking score for the compound hits from the identification nodes and pathway mapping nodes.

The mzLogic algorithm can provide a ranking score for the various database search results when an unknown compound has available data-dependent MS2 scans and similarity results from an mzCloud similarity search.

Note The ranking score is not a probability score. It is only a measure of how similar the fragmentation spectra for a putative compound are to closely matching spectra in the mzCloud spectral database.

During data processing, the mzLogic node does the following:

- 1. Runs a forward search and a reverse search using the mzCloud service.
- 2. For compounds that have available MS2 scans, scores all the structure candidates (or the specified maximum number of candidates) from the attached input nodes.

Note The following nodes can supply structures to the Apply mzLogic node: Search ChemSpider, Search Mass Lists, Map to BioCyc Pathways, and Map to Metabolika Pathways.

The mass lists that you select for the Search Mass Lists node must include structures. The Endogenous Metabolites database 4400 Compounds.masslist file does not include structures.

- 3. Adds the following columns to the result tables:
 - Adds the #Similarity Results column to the Compounds table. By default, this column is hidden.
 - Adds the mzLogic Score column to the following related tables as applicable: ChemSpider Results, Mass List Search Results, Metabolika Results, and BioCyc Results.

Running an mzLogic Analysis from the mzLogic Analysis View

If the analysis for a result file did not include the Apply mzLogic node, or if you suspect that the online mzCloud spectral database now includes more spectral data, use the mzLogic Analysis view to rank the putative structures for your unknown compounds.

Because the mzLogic Analysis view includes a link to the ChemSpider database where you can select putative structures, the result file does not need to include structure results.

Related Topics

• Running an mzLogic Analysis

Managing the Start Page's Recent File Lists

The right side of the Start Page displays lists of the most recently opened study and result files. You can clear the recent file lists, remove files from these lists, keep a specific file at the top of each list, or open the folder where a specific study file or result file resides.

See these topics:

- Keeping a File Name at the Top of a Recent File List
- Removing Files from the Recent File Lists
- Exploring File Folders from the Recent File Lists

Keeping a File Name at the Top of a Recent File List

To keep a study name or result file name at the top of its list on the Start page

Click the pin icon to the left of the name.

The orientation of the pin changes from \rightarrow to \uparrow .

Removing Files from the Recent File Lists

You can remove individual files from the recent file lists or completely clear each list.

- * To remove files from the recent files lists
 - To remove a single file, right-click the file and choose Remove From List.

-or-

• To clear a list, right-click any file in the list and choose Clear List.

Note For information about using the File menu commands to manage the recent file lists, see "Application Menu Bar" on page 7.

Exploring File Folders from the Recent File Lists

You can access the file folder for a specific study or result file from the recent files list.

* To explore the contents of a folder for a study file or result file

Under the recent files list, right-click a study file or a result file and choose **Explore Path**.

Windows Explorer opens the folder for the selected study file or result file.

Auto-Hiding the Start Page, the Chromatograms View, and the Mass Spectrum View

To make more space for other views, you can auto-hide the Start Page and the Chromatograms and Mass Spectrum views. The auto-hide features collapses each of these items down to a small tab.

To use the auto-hide feature

1. Right-click the page tab and choose **Dockable**.

The Auto Hide command becomes available for the Start Page and the Chromatograms and Mass Spectrum views. This command remains unavailable for all other dockable pages.

2. Right-click the page tab and choose Auto Hide.

The tab changes to a vertical tab on the left of the application window.

- 3. To view the hidden page, click its tab.
- 4. To hide the page, click anywhere in the application window outside the page borders.

Working with the Tabbed Documents

A tabbed document is a top-level tabbed page below the toolbar in the Compound Discoverer window.

The following pages are tabbed documents in the application window:

- Start Page (Start Page ×)
- Study files (*Study Name* ×)
- Result files (S Result File Name ×)
- Lists and Libraries manager (**11 Lists & Libraries** ×)
- Configuration (Sconfiguration ×)

- Job Queue (Z Job Queue ×)
- Report Templates (**S** *Report Template Name*)
- License Manager (<u>Q</u> License Manager ×)

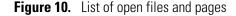
Note Excluding the Start Page and Job Queue, views that open from the View menu or toolbar are not tabbed documents—that is, when they are open, they are not listed in the open files.

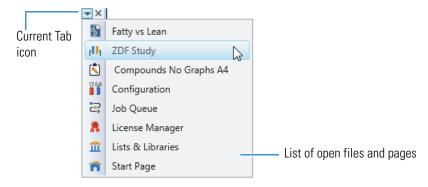
To display, dock, or hide the tabbed documents, see these topics:

- Tabbed Documents List
- Opening a Hidden Tabbed Document
- Tab Groups
- Shortcut Menu Commands for Tabbed Documents

Tabbed Documents List

Although you can have all the tabbed documents open simultaneously, including multiple study files and result files, the number of tabs that the application can display is limited by the monitor size. As you open more files than the monitor can display, the tabs begin to disappear from view in the order that you opened the files. To indicate that one or more tabs are hidden, the Current Tabs icon changes from \checkmark to \checkmark (Figure 10).





Opening a Hidden Tabbed Document

* To display an open tabbed document when its tab is hidden

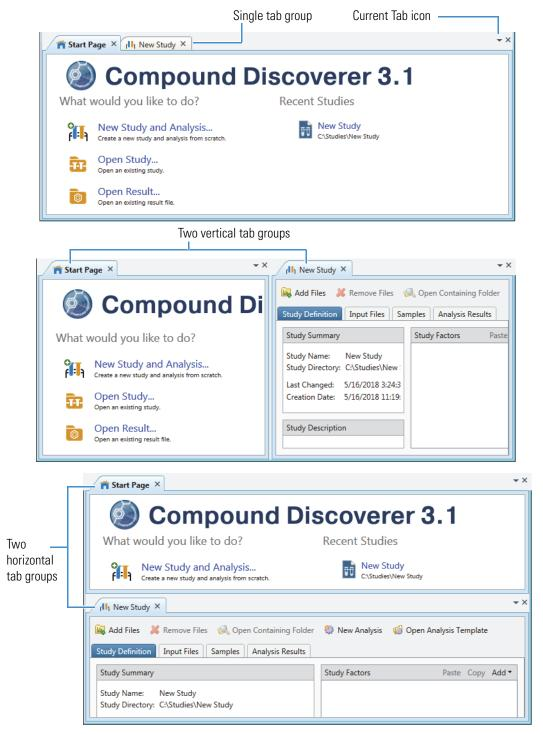
Click the **Current Tabs** icon, $\mathbf{\nabla}$, to display a list of open files. Then, select the appropriate tabbed page from the list.

The selected page becomes active.

Tab Groups

When you have two or more tabbed documents open in the same tab group, you can create more tab groups. Each tab group has its own Current Tabs icon, \checkmark . Figure 11 shows tab group examples.

Figure 11. Orientation of tab groups



Shortcut Menu Commands for Tabbed Documents

Table 7 describes the shortcut menu commands that control the tabbed document properties.

Table 7. Shortcut menu for tabbed documents

Command	Description	
Dockable	Available for the Start Page and the result file views. This command is not available for the Job Queue page.	
	Activates the Auto Hide command.	
Tabbed Document	Makes the page a tabbed document.	
	Available for the Start Page, Job Queue page, License Manager page, library pages, study pages, result file views, and report template pages.	
Auto Hide	Available for the Start Page and the views in the View menu wher these pages are dockable windows. This command is not available for the Job Queue page.	
	Hides the page while leaving the tab visible. Clicking the tab opens the page. Clicking outside the page closes the page if more than one tabbed document is open. The location of the tab depends on the tabbed document's position in the application window.	
Hide	Closes the tabbed document.	
Move to Previous Tab	Changes the position of the tabbed document.	
Group	Available only when the application window contains two or more tabbed groups.	
Move to Next Tab	Changes the position of the tabbed document.	
Group	This command is available only when the application window contains two or more tabbed groups.	
New Horizontal Tab Group	Moves the selected tabbed document to a new horizontal tab group.	
	Available only when there are two or more tabbed documents tha belong to the same tab group in the application window. Each tab group has its own Current Tab icon.	
New Vertical Tab	Moves the selected tabbed document to a new vertical tab group.	
Group	Available only when there are two or more tabbed documents tha belong to the same tab group in the application window. Each tal group has its own Current Tab icon.	

Rearranging the Tabbed Documents and Graphical Views

To make the best use of your screen space, rearrange the graphical views and tabbed documents as follows:

- Use the mouse pointer to move the graphical views that are available when you open a result file from one dock position to another dock position or to a second monitor.
- Use the mouse pointer or the shortcut menu commands to rearrange the tabbed documents within the application window.

To rearrange the tabbed documents and result page views, follow these topics as needed:

- Moving a Graphical View to Another Monitor
- Using the Mouse Pointer to Rearrange Tabbed Documents and Graphical Views
- Using the Shortcut Menu Commands to Rearrange the Tabbed Documents

Moving a Graphical View to Another Monitor

The graphical views that are available for a specific result file depend on the processing workflow used to generate the file.

* To move a graphical view to another monitor

Use the mouse pointer to drag the view by its title bar.

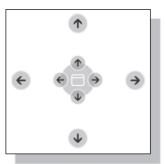
Using the Mouse Pointer to Rearrange Tabbed Documents and Graphical Views

✤ To move a graphical view or a tabbed document

1. Drag the view by its title bar or the document by its tab.

As you drag a view or a tabbed page, a guide tool appears. The guide tool consists of four directional arrows (inner arrows) that are arranged in a diamond pattern around a central circle. In addition to the guide tool, a directional arrow (outer arrows) appears in the middle of each of the window's four edges.

Figure 12. Guide tool



2. To dock a view or tabbed page in a new position, align the pointer with the appropriate directional arrow, and then release the mouse button.

Table 8. Rearranging the views and tabbed pages

To do this	Do the following
Move the selected view above the second view.	Drag the pointer to the inner up arrow, <u>(</u> .
Move the tabbed page to a horizontal group above the current group.	
Move the selected view below the second view.	Drag the pointer to the inner down arrow, 😱.
Move the tabbed page to a horizontal group below the current group.	
Move the selected view to the left of the second view.	Drag the pointer to the inner left arrow,).
Move the tabbed page to a group on the left.	
Move the selected view to the right of the second view.	Drag the pointer to inner right arrow, 💽 .
Move the tabbed page to a group on the right.	
Make both views tabbed.	Drag the title bar to the tabs icon.
	The application displays the first view and creates a tab for the second view.
Move the selected view to the top or bottom of the window.	Drag the pointer to the outer top arrow or the outer bottom arrow.
Move the tabbed page to the top or bottom of the window.	
Move the first view to the left side or right side of the window.	Drag the pointer to the outer arrow at the left of the window or the outer arrow at the right of the
Move the tabbed page to the left side or right side of the window.	window.

Using the Shortcut Menu Commands to Rearrange the Tabbed Documents

✤ To move a tabbed document to another group

Right-click the document and choose one of these commands:

- Move to Next Tab Group
- Move to Previous Tab Group
- New Horizontal Tab Group
- New Vertical Tab Group

Related Topics

Shortcut Menu Commands for Tabbed Documents

1 Introduction

Rearranging the Tabbed Documents and Graphical Views

2

Getting Started

These topics summarize the application setup, data analysis, data review, and reporting processes for the Compound Discoverer application. To get started with the Compound Discoverer application, follow these topics in order.

Contents

- 1. Setting Up the Lists and Libraries and Checking Access to the Online Databases
- 2. Setting Up a Study and an Analysis and Processing a Data Set
- 3. Opening the Result File, Reviewing the Data, and Creating a Report

Setting Up the Lists and Libraries and Checking Access to the Online Databases

To set up your lists and libraries and test your computer's access to the online databases, follow these topics as needed.

- Setting Up the Expected Compounds Library for a Targeted Analyses
- Checking the Access to the Online Databases for Identifying Unknown Compounds

Note While the application comes with a local mzVault library and several mass lists, you can also import your own custom spectral libraries and mass lists.

Related Topics

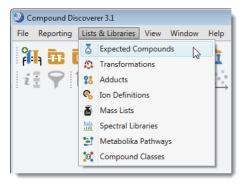
- "Using the Lists and Libraries Manager" on page 497
- "Testing Communication to the Online Databases" on page 611

Setting Up the Expected Compounds Library for a Targeted Analyses

To identify the reaction products for specific parent compounds, you must add these parent compounds to the Expected Compounds library.

- * To add a compound to the Expected Compounds library
- 1. From the menu bar, choose Lists & Libraries > Expected Compounds.

Figure 13. Lists & Libraries menu



2. Click New.

The Compound Editor opens.

Figure 14. Compound Editor

② Compound Editor	
N 🖓 🖓 🗸 🗸 🖉 🖌 🖉	<u> </u> () (Ф (
	_
Name:	
Name:	
Description:	
Elemental composition:	
Molecular weight:	
0.00000	
ChemSpider	Cancel

3. At the bottom of the editor, click **ChemSpider**.

The ChemSpider Search window opens.

4. Enter the name of a compound and click **Search**.

Figure 15. ChemSpider Search window with a result for caffeine

CSID Name Formula Molecular Weight # References Structure = * Aa * = * = * Aa *
= • Aa • = • = • Aa •
2424 Caffeine C8 H10 N4 O2 194.0803755 13026

5. Click Select.

The ChemSpider Search window closes.

6. In the Compound Editor, click Save.

The new compound appears in the library.

Figure 16. Expected Compounds library with caffeine (new entry) and omeprazole (existing entry)

Ocmpound Discoverer 3.	1.0					
<u>File R</u> eporting Lists & L	ibraries <u>V</u> ie	ew <u>W</u> indow	<u>H</u> elp			
îi 🖬 🛅 🖥						
i E 🌱 🕰 🖞	L, C,			V, Ľ (0. 😂 🖻	백 백 장 중 🖨 🗄 🕄
😭 Start Page 🗙 🏦 L	ists & Libra	ries ×				- ×
5	New		Edit	Delete	Import	xport All
Expected Compounds	₽	Name 🔺	Description 🔺	Elemental Composition 🔺	Molecular Weight [Da] 🔺	Structure
82		<u>A</u> a •	<u>A</u> a +	<u>A</u> a •		<u>A</u> a •
Transformations	1 🗢	caffeine		C8 H10 N4 O2	194.08038	
28						O N N N
Adducts	2 🖙	omeprazole		C17 H19 N3 O3 S	345.11471	
Ion Definitions						~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~
	4				Ш	×

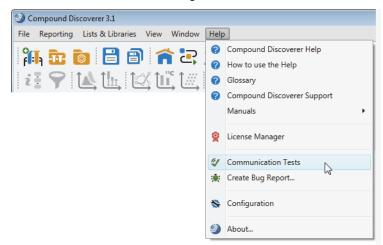
Related Topics

• "Modifying the Expected Compounds List" on page 498

Checking the Access to the Online Databases for Identifying Unknown Compounds

Check the computer's access to the KEGG Pathways, BioCyc Pathways, mzCloud, and ChemSpider databases.

- * To check the computer's access to the online databases
- 1. From the menu bar, choose **Help > Communication Tests**.



- 2. Click the database's tab.
- 3. Click Run Tests.

Figure 17. mzCloud communication tests

🕼 Communication Tests	×
BioCyc KEGGPathway mzCloud ChemSpider	
Checking service availability Succeeded	
Checking credentials Succeeded	
Succeeded	
Succeeded	
Checking batch search Succeeded	
6/22/2019 4:24:13 PM mzCloud: Checking batch search - Starting batch search using generated test spectrum	
-> Succeeded	
Clear Results Copy Info Run Tests Clos	e

4. (Optional) To use the BioCyc database, set up a BioCyc subscription.

Related Topics

- "Running the Communication Tests" on page 612
- "Setting Up a BioCyc Account or Subscription" on page 575

Setting Up a Study and an Analysis and Processing a Data Set

A study is where you define your samples. An analysis is where you specify how you want to analyze the raw data (with a selected processing workflow), the samples (raw data files) that you want to process, (optional) the sample groups and ratios for a differential analysis, and the name of the result file for storing the processed data.

Follow these topics in order:

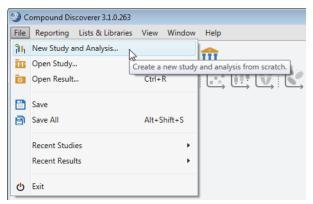
- 1. Setting Up a New Study and a New Analysis by Using the Wizard
- 2. Troubleshooting an Analysis
- 3. Submitting a Run to the Job Queue

Setting Up a New Study and a New Analysis by Using the Wizard

To set up an new study and a new analysis

1. From the menu bar, choose File > New Study and Analysis.

Figure 18. File menu



The New Study and Analysis Wizard opens. To display instructions for each page of the wizard, click the **Show Description** icon, **?**.

- 2. On the Study Name and Processing Workflow page, do the following:
 - a. In the Study Name box, type a name for the study.

Note The application creates a study file and a study folder with this name. The application stores the study file and all the result files that you generate by running analyses within the study in the study folder.

b. Click the browse icon to the right of the Studies Folder box, and then browse to the folder where you want to store your studies.

Note You can store all your studies in this top-level folder—that is, you do not need to create a new top-level Studies folder each time you create a new study.

When you create a study, the application creates a study folder with the same name as the study and stores this folder in the top-level Studies folder.

- c. In the Processing area, open the Workflow list, and then select a processing workflow template for your area of study:
 - Degradant ID

• Environmental

- Forensics
- Natural Products

- Extractables & Leachables
- Metabolism (MetID)
- Polymers
- Metabolomics
- Stable Isotope Labeling
- d. Read the description of the processing workflow.

Figure 19 shows the selection of a processing workflow that identifies detected compounds by searching online databases.

tudy Name and Director	y Structure	
Study Name:	Caffeine Study	
Studies Folder:	C:\Studies	
Study Template File:	(Optional)	
Description:	(Optional)	
Processing		
Workflow:	WorkflowTemplates \ Metabolomics \ Untargeted Metabolomics using Online Databases, mzLogic, and Molecular Networks 💌	
Workflow Description:	Untargeted Metabolomics workflow: Find and identify the differences between samples. - Performs retention time alignment, unknown compound detection, and compound grouping across all samples. Predicts elemental compositions for all compounds, fills gaps across all samples, and hides chemical background (using Blank	

- e. Click Next.
- 3. On the Input File Selection page, do the following:
 - a. Click Add Files.

The Add Files dialog box opens.

- b. Browse to and select the raw data files that you want to add to the study, and then click **Open**.
- c. Click Next.

Tip To quickly learn the basic workflow for the application—from creating a new study to reviewing a result file—skip step 4 and step 5, as setting up the study variables can be time consuming. You can set up the study factors and the sample groups for your sample set after you become more familiar with the application.

For a guided tutorial to setting up a study, including setting up the study factors for a sample set, follow the tutorial for your area of study. You can access these tutorials by choosing Help > Manuals from the application menu bar.

4. (Optional) On the Input File Characterization page, define and assign the study factors for the sample set and assign the sample types. Then, click **Next**.

Tip To define the study factors and characterize the input files, follow the instructions in the left pane or press the F1 key, **1**, on your computer keyboard to open the Help to the **Characterizing New Input Files** topic.

5. (Optional) On the Sample Groups and Ratios page, set up the sample groups and ratios for a differential analysis. Then, click **Next**.

Tip To set up the sample groups and ratios, follow the instructions in the left pane or click anywhere on the Sample Groups and Ratios page, and then press the F1 key, , on your computer keyboard to open the Help to the **Setting Up the Sample Groups and Ratios for a New Analysis** topic.

6. Click **Finish** to exit the wizard.

The study opens as a tabbed document at the left of the application window, and the Analysis pane opens to the right of the study.

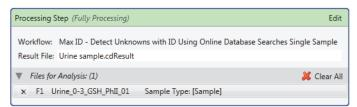
The study consists of four tabbed pages: Study Definition, Input Files, Samples, and Analysis Results. The two tabs (Grouping & Ratios and Workflows) to the right of the study pages are part of the analysis—that is, if you close the Analysis pane, these tabs also close.

The Analysis pane lists the selected processing workflow, the name of the result file, and the selected raw data files. If the analysis is valid, the Run button is green. Otherwise, the Run button is unavailable.

		Stu	udy pages		_	Analy	vsis pages	5	ļ	Analysis pane		Run bı	itton	
2	ompour	nd Discoverer 3.1												×
<u>F</u> ile	<u>R</u> epor	ting <u>L</u> ists & Librarie	es <u>V</u> iew <u>W</u> indo	ow <u>H</u> elp										
f f	4 0 2 5) 🟫 🗉				Zé	1	<u>)</u> 124		· & [₽ (1)	(
1		Page × II Caffei		-	ŕ									. ×
	Add Fil	es 💥 Remove File	s 🔍 Open Co	ntaining Folde	L	New Analysis	🕼 Open		is Template Analysis		As Bate	th 💰 Run	E Save	×
	r ID 4		Jampies And	File Type	U	Sample Informat		<u> </u>	Analysis		AS DOLL	n 🤤 Kun	Jave	
		•		21	•		•		Processing	Step (Fully Processing)			E	dit
	F1	Urine_0-3_GSH_PhI	[_01	.raw		Sample Type: [Sa	ample]		Workflow	 Untargeted Metabolom mzLogic, and Molecular Urine 0-3 GSH_PhII 01. 	Networks	e Database	s,	
											cakesuit			
										or Analysis: (1)			样 Clear	All
									× F1	Urine_0-3_GSH_PhII_01	Sample Type	: [Sample]		
	Show D	etails												
											th	efault file e result f :dResult)		or

Figure 20. Study page and Analysis pane

7. (Optional) In the Result File box of the Analysis pane, rename the result file.



Note By default, the application uses the name of the first input file as the result file name.

Related Topics

• "Starting a New Study and Setting Up an Analysis" on page 67

Troubleshooting an Analysis

* To troubleshoot the analysis when the Run button is unavailable

1. In the Analysis pane, point to the Caution symbol, 🔔.

A pop-up message lists the missing analysis items.

If you are working with one of the defined processing workflows, the most common errors are not selecting the following items for the workflow nodes:

- Compounds for the Generate Expected Compounds and Create FISh Trace nodes
- A mass list for the Search Mass Lists node
- A spectral database for the Search mzVault node
- An isotope pattern for the Pattern Scoring node
- The isotope ratios for the Create Pattern Trace node
- A Compound Class fragment list for the Compound Class Scoring node
- 2. Click the Workflows tab and fix the analysis errors until the Caution symbol disappears.

To fix a missing parameter error for a node (**9**), select the node in the Workflow Tree pane on the Workflows page. Then, on the Parameters page for the node, select the missing item.

For example, if you have not specified any compounds for the Generate Expected Compounds node, select the node in the Workflow Tree area, and then select one or more compounds from the Compound list.

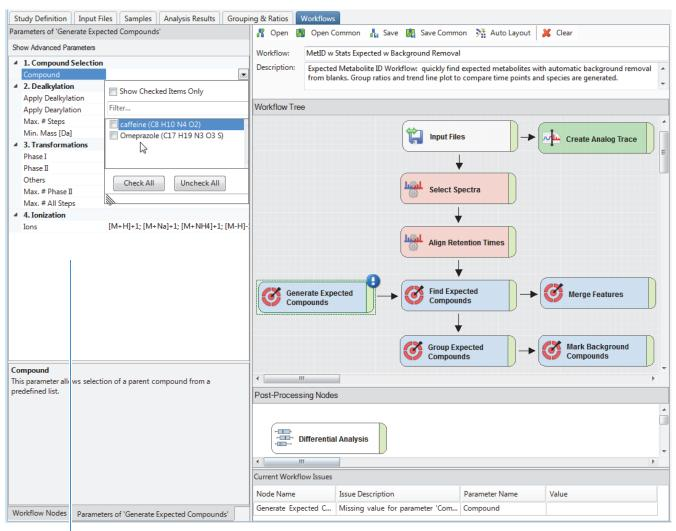
Note The Current Workflow Issues table at the bottom of the Workflows page also lists all the processing workflow's issues. This table does not list other validation issues, such as no input files in the Files for Analysis area of the Analysis pane.

Tip In Compound Discoverer 3.1 or later application, you can leave the analysis open as you update any of the lists under Lists & Libraries.

Figure 21 shows a processing workflow template for a targeted analysis. For this template, you must select one or more compounds for the Generate Expected Compounds node.

2 Getting Started

Setting Up a Study and an Analysis and Processing a Data Set





 Parameters for the Generate Expected Compounds node

Figure 22 shows the Max ID template for Metabolomics studies. This template is only suitable for a single sample, as it does not include the Align Retention Times node, which is required for aligning the detected compounds across all the input files. Because the template is designed to process a single sample, it does not include the Differential Analysis node under Post Processing Nodes.

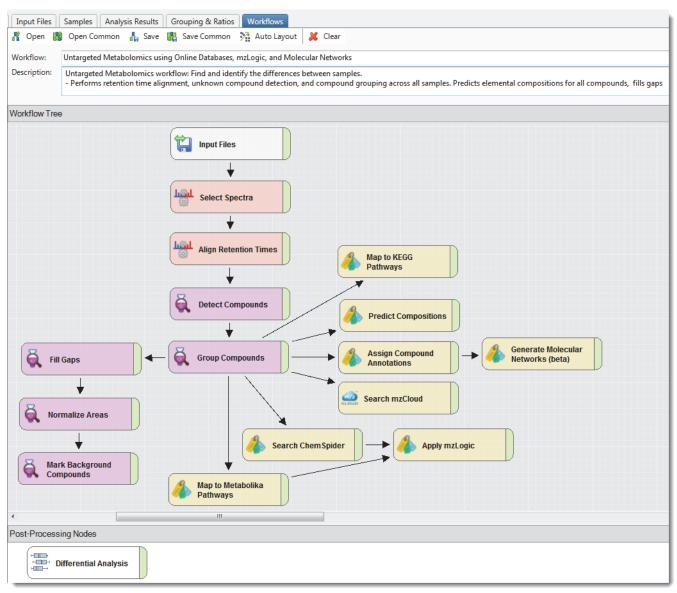
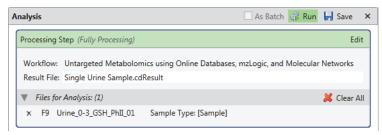


Figure 22. Untargeted processing workflow with mzLogic and Molecular Networks

Submitting a Run to the Job Queue

To submit a run to the Job Queue

- 1. In the Analysis pane, do the following:
 - a. Rename the result file as appropriate.



Note By default, the application gives the result file the same name as the first input file.

- b. Click Run.
- 2. If the Analysis Validation Issue prompt opens, decide whether to ignore the issue or cancel the run.

Tip Not defining the ratios on the Grouping & Ratios page when the processing workflow includes the Differential Analysis node is one of the most common validation issues. If you do not want the processing workflow to perform a differential analysis, click **Ignore**. Otherwise, click **Cancel**, set up the sample groups and ratios, and then click **Run** again.

If there are no validation issues or you click Ignore to continue submitting the job, the Job Queue page opens.

Opening the Result File, Reviewing the Data, and Creating a Report

Note When a graphical view or result table is the active item, pressing the F1 key opens the Help topic for the specific view or table.

To review the processed data and create a report, follow these topics in order:

- 1. Opening the Result File from the Job Queue
- 2. Reviewing the Compounds that the Analysis Detected or Found
- 3. Displaying Only the Items of Interest in a Result Table
- 4. To create a report, follow either of these topics:
 - Exporting a List of Compounds to a Spreadsheet
 - Printing a Report Using a Defined Template

Opening the Result File from the Job Queue

* To open a result file

- 1. If the Job Queue page is closed, choose **View > Job Queue** from the application toolbar.
- 2. On the Job Queue page, double-click the completed job.

The result file opens as a tabbed page.

Reviewing the Compounds that the Analysis Detected or Found

✤ To review the processed data

1. Select a compound of interest in the Compounds table or the Expected Compounds table.

The respective XIC traces appear in the Chromatograms view and the best MS1 spectrum appears in the Mass Spectrum view.

Figure 23 shows the result page for a single sample that was processed with the following workflow template: Untargeted Metabolomics using Online Databases, mzLogic, and Molecular Networks.

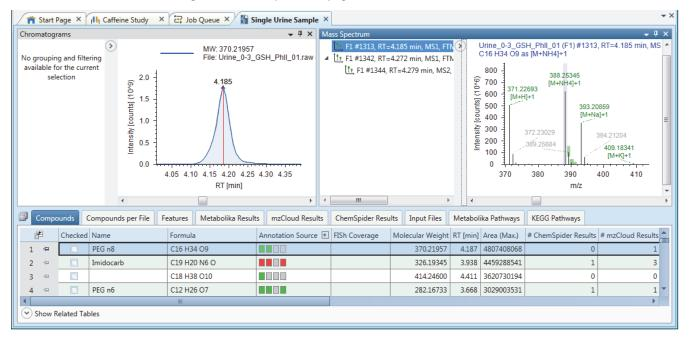


Figure 23. Example result page

2. Review the various result tables and open other views as applicable.

Related Topics

• "Reviewing the Analysis Results" on page 247

Displaying Only the Items of Interest in a Result Table

Use the Result Filters view to display only the compounds of interest in the Compounds table (for an untargeted analysis) or the Expected Compounds table (for a targeted analysis).

* To display only the checked compounds in the table

- 1. Open a result file and click the **Compounds** tab or the **Expected Compounds** tab.
- 2. In the result table that you selected, select the compounds of interest by selecting their respective check boxes in the Checked column.
- 3. From the menu bar, choose **View > Result Filters**.

Figure 24. View menu

File F	Report	ing Lis	ts & Libra	aries	View	Window Help	
: ក្រ	н- e	: 🗎	🖻 : 👩	3	n	Start Page	Ctrl+H (
1	Start F	age ×	rll ₁ Caf	feine S	2	Job Queue	Ctrl+J
Chron	natogr	ams			i	Result Summary	ec
			>)	9	Result Filters	1
	able fo	g and filt or the cu action		1 (1 0^9)	[<u>▲</u> [<u>⊥</u> ,	Chromatograms Mass Spectrum	1
				Intensity (counts) (10^0)	<mark>⊠</mark> ≣	Trend Chart Isotopologues Distribution Chart Mass Defect Plot	
				*	[<u>.</u> [<u>.</u> [⊻	Result Charts Descriptive Statistics Differential Analysis	
	Compo	ounds	Compou	nds p	2	Principal Component Analysis	mS
f	ŧ	Checke	d Name		\mathbb{Z}	Partial Least Squares - Discriminant Analysis	Co
1	Þ		PEG n8	1	Ľ	Hierarchical Cluster Analysis	
2 3	ф Ф		Imidoc	arb	(Ö)	Retention Time Corrections Compound Area Corrections	
4	÷		PEG n6		<u>_</u>	Compound Area Corrections	
5	12		E IL I		Þ.	Metabolika Pathways	-
6 7	4 4		5-Hydr	oxyor	Ū,	BioCyc Pathways	-
8	Ę.				¢,	KEGG Pathways	
9	- -	1	Caffein	e	٢	mzLogic Analysis	
						FISh Scoring Queue	
⊘s	how R	elated Ta	ables		-		

The Result Filters view opens as a floating window.

4. In the Result Filters view (Figure 25), under Compounds in the right pane, click **Add Property**, and then select **Checked** from the drop-down properties list.

② Result Filters				
ON Compounds ON Compounds per File ON Features ON Matabolika Results ON Metabolika Results ON ChemSpider Results ON Input Files ON Metabolika Pathways ON Metabolika Pathways	AN OR # 4 # C # K # M # n # s An Are Are Baa Gre For Gra Gra	Add aroup property) D dducts hemSpider Results EGG Pathways EGG Pathways testabolika Pathways nzCloud Results miliarity Results molarity Results nocurce a a (Max.) kground scked h Coverage mula p Status sup Areas SG Pathways		II Clear Apply Filters
	Me Mc MS	tabolika Pathways Iecular Weight	Ŧ	

Figure 25. Adding the Checked property for filtering the Compounds table

- 5. To apply the filter to the Compounds table, click **Apply Filters**.
- 6. In the Compounds table, only the checked compounds remain visible.

		Compo	unds 🏆	Compounds per File	- Features	Metabolika	Results	mzCloud Resu	lts
[É	₽	Checked	Name	Formula		Annotati	ion Source	+
	1	- 1 2	V	5-Hydroxyomeprazole	C17 H19 N3 C	04 S		l	
	2	÷	V	Caffeine	C8 H10 N4 O2	2			

Exporting a List of Compounds to a Spreadsheet

- To export the compounds data to a spreadsheet
- 1. Use the Result Filters view to filter the data or select the check boxes for the compounds of interest.
- 2. Right-click the Compounds table or the Expected Compounds table and choose the appropriate format:
 - a. To export all the displayed columns, choose **Export > As Excel**.

Copy With Headers	Ctrl+C	ľ	
Сору			
Clear Selection		L	
Cell Selection Mode		L	
Enable Column Fixing		l	
Collapse All Column Headers			
Expand All Column Headers			
Check Selected	•		
Check All	•		
Uncheck Selected	•		
Uncheck All	•		
Remove All Checkmarks in All Tables			
Edit Compound Annotation			
Clear Compound Annotation		L	
Apply FISh Scoring			
Molecular Networks (beta)		L	
Molecular Networks (Deta)		L	
Export	•		As Plain Text
			As Excel
			As Xcalibur Inclusion/Exclusion List
			As TraceFinder List
			As mzVault Library
			Add Compound to Existing mzVau
			As Mass List
			Add Selected Compounds to Existi

Figure 26. Shortcut menu for the Compounds table

The Export to Excel dialog box opens.

- b. Select the result tables of interest.
- c. To export only the checked compounds, select the Checked Items Only check box.

C:\Users\Public\Documents\Single Urine Sample.xlsx		
Items and related tables to be exported Level 1: Compounds Level 2: Level 3:	▼ ▼ ▼	Options Checked items only Open file after export

- d. In the Path box, browse to the folder where you want to store the spreadsheet.
- e. Click **Export**.

The application exports the spreadsheet to the selected folder.

Printing a Report Using a Defined Template

* To print a report using a defined report template

- 1. Open the Expected Compounds table or the Compounds table.
- 2. Filter the table to display only the compounds of interest.
- 3. From the application menu bar, choose **Reporting > Create Report**.
- 4. In the Open Report Design Template dialog box, select one of the Compounds report templates or one of the Expected Compounds templates.

Figure 27. Report templates

Irganize ▼ New folder					•	
Name	Date modified	Туре	Size			
Compounds No Graphs A4.cdReportTemplate	7/4/2019 9:55 AM	CDREPORTTEMPLATE File	28 KB			
Compounds with Graphs A4.cdReportTemplate	7/4/2019 9:56 AM	CDREPORTTEMPLATE File	45 KB			
Expected Compounds No Graphs A4.cdReportTemplate	7/4/2019 9:57 AM	CDREPORTTEMPLATE File	30 KB			
Expected Compounds per File No Graphs A4.cdReportTemplate	7/4/2019 9:57 AM	CDREPORTTEMPLATE File	31 KB			
Expected Compounds per File with Graphs A4.cdReportTemplate	7/4/2019 9:58 AM	CDREPORTTEMPLATE File	48 KB			
Expected Compounds with Graphs A4.cdReportTemplate	7/4/2019 9:59 AM	CDREPORTTEMPLATE File	47 KB			
Expected Compounds with Structures No Graphs A4.cdReportTemplate	7/4/2019 10:00 AM	CDREPORTTEMPLATE File	30 KB			
Expected Compounds with Structures with Graphs A4.cdReportTemplate	7/4/2019 10:01 AM	CDREPORTTEMPLATE File	47 KB			
File name: Compounds with Graphs A4.cdReportTemplate			-	CD Report Te	mplate Fi	le (*.cdR

5. Click **Open**.

The report resolution page opens and the application resolves the data with the template (Figure 28).

Opening the Result File, Reviewing the Data, and Creating a Report

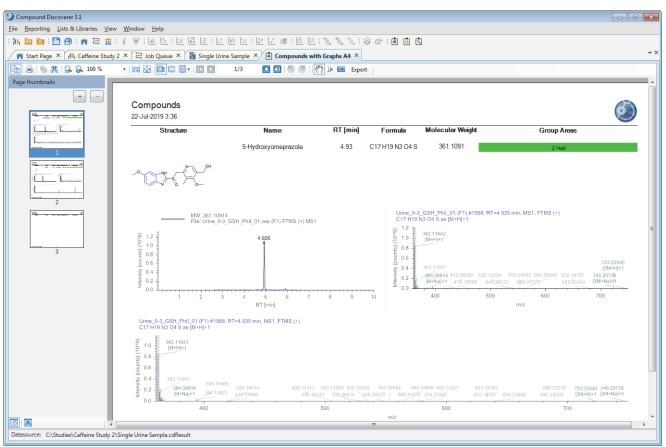


Figure 28. Resolved report

- 6. On the report resolution page, click the **Print** icon, (a), in the toolbar to print the report.
- 7. In the Print dialog box, do the following:
 - a. Select the appropriate printer and the page range that you want to print.

Note The report templates that come with the application default to printing on A4 paper.

- b. If you are not printing on A4 paper, change the printer setting.
- c. Click **OK** to print the report.

Related Topics

• "Generating a Report with an Existing Report Template" on page 332

3

Starting a New Study and Setting Up an Analysis

To create a new study and optionally to set up an analysis, see these topics.

Contents

- About Studies and Analyses
- Data Storage Architecture
- Starting the New Study and Analysis Wizard
- Setting Up the Study Folders and Selecting a Processing Workflow
- Adding Input Files to a New Study
- Characterizing New Input Files
- Using the Regular Expression Builder to Extract the Study Factors
- Setting Up the Sample Groups and Ratios for a New Analysis
- Preparing to Submit a Run to the Job Queue
- About Accessing the Wizard Help

About Studies and Analyses

In the Compound Discoverer application, you start analyses (data processing) from within the study environment.

For information about studies and analyses, see these topics:

- Studies
- Input Files
- Study Factors
- Sample Types
- Analyses

Studies

A study includes a list of input files with their location, the sample information for each file, and a list of the analyses run within the study. The sample information includes the file name, the study factor values, and the sample type for each sample. Creating a new study automatically creates a study file (.cdStudy) for storing the study information and a study folder for storing the study file and any result files (.cdResult) generated by running analyses within the study. The study file and its folder share the same name (see "Data Storage Architecture" on page 70). **Input Files** Input files are the raw data files acquired from an LC/MS/MS experiment. The file format for raw data files acquired with a Thermo Scientific MS is RAW. To predict the elemental compositions of the sample components, the raw data must include high-resolution accurate-mass MS1 scans. To search the mass spectral libraries for matching fragmentation spectra, the raw data must include data-dependent MS2 or higher scans. **Study Factors** A study factor is an experimental variable that might have a statistically significant effect on the sample population being studied. There are three types of study factors: Categorical Factors include non-quantifiable categories such as organism, matrix, tissue, gender, and so on. • Numerical Factors include quantifiable variables such as time, amount, concentration, and so on. You can assign a unit to a numeric study factor; however, the unit is only a label and is not used in any calculations. • Biological Replicate Factors include non-quantifiable categories with samples from different entities of the same type under the same conditions and provide a measure of the variability associated with these conditions. Use the biological replicate factor for nested experiments. See Appendix A, "Experiment Design for Comparison Statistics." You can add only one biological replicate factor to a study. **Sample Types**

Table 9 describes the sample types that the application supports and how the it uses these sample types.

Sample type	Application use
Sample ^a	Detects the compounds in the sample.
Control ^a	Detects the compounds in the sample.
Blank	Detects the compounds in the sample. When the processing workflow includes one or both Mark Background Compounds nodes, marks these components as background compounds.
Identification Only	Does not report the chromatographic peak areas for the sample's compounds. Uses the sample's fragmentation scans for component identification when the processing workflow includes the Group Compounds node.
Quality Control	Detects the compounds in the sample. Pools the QC samples to determine a group area for each detected compound for area normalization.
Standard ^a	Detects the compounds in the sample.
Labeled	Determines the formulas for the compounds in the labeled samples.

^a The application attributes the same functionality to the Sample, Control, and Standard sample types. You can use the Control and Standard sample types to label your control or standard samples; that is, you can use these sample types as an additional study variable for grouping.

Related Topics

• Selecting the Sample Types

Analyses

An analysis includes the processing workflow that extracts information from the raw data for each sample, such as the chromatographic peak areas and mass spectra for known and unknown compounds. It also includes the group ratios required to run a differential analysis, a list of input files, and the name of the result file where the processing results reside.

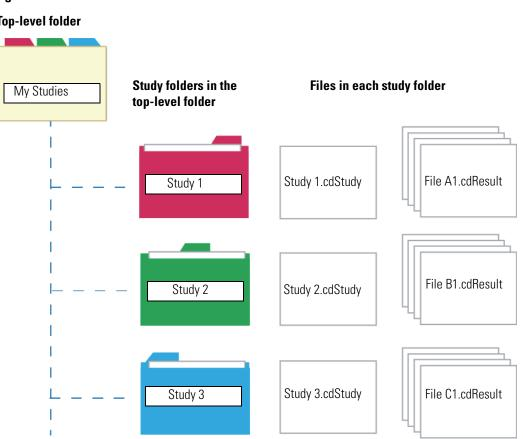
Related Topics

- Creating, Running, and Reprocessing Analyses
- Understanding Processing Workflows
- Reviewing the Analysis Results

Data Storage Architecture

Figure 29 shows the hierarchy of the application folders. The studies folder is the top-level folder for all your studies or a particular set of studies. Each study folder within the studies folder holds one study file (.cdStudy) and one or more result files (.cdResult).

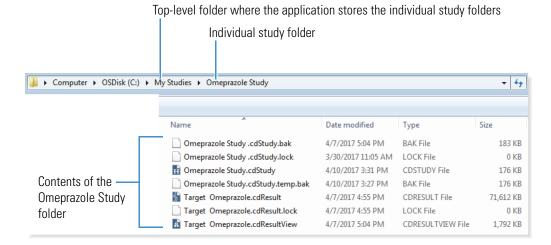
Note You can create more than one top-level folder for your studies. Each time you open the New Study and Analysis Wizard, it opens to the last opened top-level folder.





Top-level folder

This figure shows an example of the directory hierarchy for a study named Omeprazole Study with one result file. The individual study folder and the study file share the same name.



Starting the New Study and Analysis Wizard

You can access the wizard from the Start Page, the File menu, or the application toolbar.

* To start the New Study and Analysis Wizard

Do one of the following:

- From the menu bar, choose File > New Study and Analysis.
- In the toolbar, click the Create a New Study and Analysis icon, 1.

-or-

• On the Start page under What Would You Like to Do?, click **New Study and Analysis**.

The New Study and Analysis wizard opens. Unlike other areas of the application, the wizard has embedded "How To" instructions.

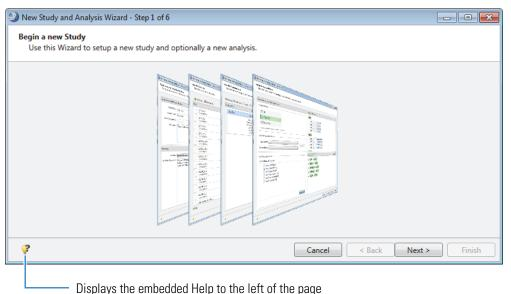


Figure 30. First page of the New Study and Analysis Wizard

Setting Up the Study Folders and Selecting a Processing Workflow

Use the Study Name and Processing Workflow page (step 2 of 6) of the New Study and Analysis Wizard to do the following:

- Name the current study and its storage folder.
- Select the top-level folder for storing all the study folders or a subset of the study folders.
- (Optional) Select a template file for the study.

Note A template file includes defined study factors.

• (Optional) Select a processing workflow for the analysis.

Note You can select a processing workflow after you create the study.

Follow these topics as needed:

- Naming a Study and Its Folder
- Creating or Selecting the Top-Level Folder for the Study
- Selecting a Study Template
- Selecting a Processing Workflow for the Analysis

After you complete this page of the wizard, go to the next page where you select the input files for the analysis.

Figure 31 shows page 2 of the wizard.

Figure 31. Study Name and Processing Workflow page of the wizard

	or this study and its folder, select the studies folder for storing all of your study cessing workflow for the current analysis.
Study Name and Directo	ny Structure
Study Name:	Lean versus Fatty
Studies Folder:	C:\Compound Discoverer Studies
Study Template File:	(Optional)
Description:	(Optional)
Processing	
Workflow:	(empty workflow)

Naming a Study and Its Folder

The application uses the specified name for the study file (.cdStudy) and the study folder where it stores the study file. In addition, the application stores the result files from analyses run within the study in the same study folder.

To name the study and its folder

On page 2 of the wizard, in the Study Name box, type a name for the new study.

For more information, see "Data Storage Architecture" on page 70.

Creating or Selecting the Top-Level Folder for the Study

The studies folder is where the application stores the individual study folders.

* To specify the top-level folder where you want to store the new study folder

On page 2 of the wizard, do the following:

• Use the current folder name and location that is displayed in the Studies Folder box.

-or-

• Browse to or create a new folder.

The new folder name and location appear in the Studies Folder box.

Selecting a Study Template

A study template is a study that includes defined study factors.

- To extract the study factors for the current study from an existing study
- 1. On page 2 of the wizard, click the browse icon, ..., next to the Study Template File box.
- 2. In the Select Template Study File dialog box, locate a study file (.cdStudy) with the appropriate study factors, and click **Open**.

The name and location of the study file appear in the Study Template File box.

Selecting a Processing Workflow for the Analysis

You can select a processing workflow on page 2 of the wizard or on the Workflows page of an analysis.

* To select a processing workflow for the current analysis

Do either of the following:

• Select a processing workflow from the Workflow list.

The Workflow list displays the processing workflow files in the Common Templates folder.

-or-

• Select a processing workflow or result file from another folder by clicking the browse icon,, next to the Workflow list.

When you select a processing workflow file, the file name appears in the Workflow list. When you select a result file, the following text appears in the Workflow list:

Imported from: File name

where File name is the file name of the result file

If the processing workflow includes a description, the description appears in the Workflow Description box below the Workflow list.

Adding Input Files to a New Study

Use the Input File Selection page (step 3 of 6) of the New Study and Analysis Wizard to select all the input files for the new study or only those input files that you want to process with the current analysis. If you click Finish before you add files to the study, the wizard saves the named study to the named folder and closes.

Note The example Xcalibur RAW files for the tutorials are on the Compound Discoverer USB key in the software media kit or in the software media that you downloaded from Flexera. The software installation process does not install the example files on your processing computer.

* To add or remove raw data files

- 1. On page 3 of the wizard, do the following:
 - a. To add files, click **Add Files**. Then, browse to the appropriate folder, select the Xcalibur RAW files of interest, and click **Open**.

Figure 32. Input File Selection page of the wizard

New Study and A	Analysis Wizard - Step 3 of 6			_	
Input File Selecti Select the inpu	ion ut files for this analysis.				
🛺 Add Files	💥 Remove Files				
Files					
0 files					
<i>?</i>		Cancel	< Back	Next >	Finish

The file names of the selected files appear in the Files box, the number of files that you selected appears below this box, and the Next button becomes available (Figure 33).

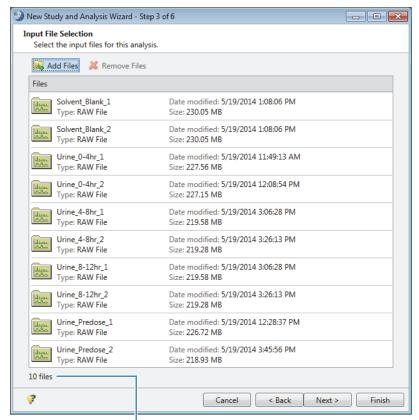


Figure 33. Input File Selection page with selected files

Number of files field

- b. To remove any of the added files, select them and click **Remove Files**.
- 2. Do one of the following:
 - Click **Next** to continue to the Input File Characterization page.
 - Click **Finish** to create the new study and close the wizard.

Characterizing New Input Files

To set up the study factors and the sample information for the study's input files, use the Input File Characterization page (step 4 of 6) of the New Study and Analysis Wizard or the Input File Characterization dialog box that opens when you add input files to an existing study.

The sample information includes the sample type and study factor values for each sample.

Note In an existing study, you can do the following:

- Use the Input File Characterization dialog box to add and edit the study factors and set up the sample information for additional input files.
- Add and edit the study factors on the Study Definition page.
- Manually select the sample types and study factor values on the Samples page.

Figure 34 shows the functional area of the Input File Characterization page or dialog box.

Figure 34. Input File Characterization page of the wizard

Opens the Extract Sample Information from Sample Names dialog box-----

Automatically assigns the defined study factor values -

Paste Copy Ac	d 🕶	Samp	ples									
Edit Y		Error	Sampl	e 🔺	File	Sample Identifier	Sample Ty	pe	Phenoty	/pe		
				•				•		*		
Lean			S1		F1	blank	Blank	*	n/a			
)			S2		F2	Pooled_ddMS2	Sample	•	n/a			
			S3		F3	ZDF_Fatty_1	Sample	*	Fatty			
				S4		F4	ZDF_Fatty_2	Sample	•	Fatty		
					S5		F5	ZDF_Fatty_3	Sample	*	Fatty	
						S6		F6	ZDF_Lean_1	Sample	•	Lean
			S7		F7	ZDF_Lean_2	Sample	•	Lean			
			S8		F8	ZDF_Lean_3	Sample	*	Lean			
	Edit × Fatty	Edit × Fatty	Edit × Fatty	Edit X Fatty Lean S1 S2 S3 S4 S5 S6 S6 S7	Edit X Fatty Lean Error Sample Error Sample S1 S2 S3 S4 S5 S6 S7	Edit × File Fatty Image: Complex of the second se	Edit X Fatty Lean S1 F1 blank S2 F3 ZDF_Fatty_1 S4 F4 ZDF_Fatty_2 S5 F6 ZDF_Fatty_3 S6 F6 ZDF_Lean_1 S7 F7	Edit X File Sample Identifier Sample Ty Fatty Image: Construction of the state of the st	Edit X Fatty Fatty Sample File Sample Identifier Sample Type B •	Edit X File Sample Identifier Sample Type Phenoty Fatty Image: Comparison of the system of the syste		

Follow these topics:

- 1. Selecting the Delimiters for Parsing the File Names
- 2. To set up the study factors, follow these topics as needed:
 - Adding Categorical Study Factors
 - Adding Numeric Study Factors

Characterizing New Input Files

- Adding Biological Replicate Study Factors
- Deleting Study Factors
- Duplicating Study Factors
- Editing Study Factors
- 3. To select or assign the study factor values to the samples (in the Samples area), follow these topics:
 - Automatically Assigning Study Factor Values
 - Manually Selecting the Study Factor Values
 - Resetting the Sample Assignments
- 4. To select the sample types, see Selecting the Sample Types.

Selecting the Delimiters for Parsing the File Names

You can select the delimiters for parsing a file name on the Input File Characterization page of the wizard or in the Input File Characterization dialog box.

- To select the delimiters that separate the factors in the file names
- 1. Select the check box or check boxes of the delimiter or delimiters for the input file names.

Delimiters: 🗹 Underscore 🔲 Hyphen 📄 Comma 📄 Space 💭 Plus 💭 Other

For example, the underscore character is the delimiter that separates the study factors from the other parts of the file name in the following file names:

Urine_0-4hr_1	Urine_4-8hr_1	Urine_8-12hr_1	Urine_Predose_1
Urine_0-4hr_2	Urine_4-8hr_2	Urine_8-12hr_2	Urine_Predose_2

In this case, the study factors are time period (0-4hr, 4-8hr, 8-12hr, or Predose) and replicate (1 or 2).

2. If the delimiter is not available, select the **Other** check box and type the delimiter character in the box.

Adding Categorical Study Factors

You can add study factors when you create a study, when you add files to an existing study, and on the Study Definition page of an existing study.

✤ To add a categorical study factor

From the menu bar at the top right of the Study Factors pane, choose Add > Categorical Factor.

The categorical factor editor appears with [new factor] automatically selected (Figure 35).

Figure 35. Categorical factor editor

AB [new factor]	Apply Cancel 🗙
Items:	
	Add Delete

- 2. Type the name of the factor, for example, **Time Period**.
- 3. For each item that you want to add to the Items list, do the following:
 - a. In the Items box (to the left of the Add button), begin typing a factor, for example,
 0-4hr.

If the file name contains a character delimiter and you selected the delimiter's check box, the editor automatically enters the appropriate text in the Items box as you start typing. Otherwise, you must type all the characters for the item.

Note The file name parsing feature is not available on the Study Definition page of an existing study, so you must type all the characters for the item.

The Add button becomes available (Figure 36).

Figure 36. Entering items in the item box

Time Period	Apply Cancel 🗙
Items:	
0-4hr	Add Delete

b. Click **Add**.

The current item appears in the Items list.

4. To save the study factor, click **Apply**.

The factor editor collapses to show only the study factor name and the Items list (Figure 37). The items appear in ascending order.

Figure 37. Named factor with a list of items

AB Time Period	Edit ×
	0-4hr 4-8hr 8-12hr Predose

Adding Numeric Study Factors

You can add study factors when you create a study, when you add files to an existing study, and on the Study Definition page of an existing study.

✤ To add a numeric factor

From the menu bar at the top right of the Study Factors pane, choose Add > Numeric Factor.

The numeric factor editor appears with [new factor] automatically selected (Figure 38).



[new factor]	Apply Cancel $ imes$
Factor Unit:	
Values:	
Add	Delete

- 2. Type a factor name to replace [new factor], for example, Replicate.
- 3. Point to the right of Factor Unit and, in the box that appears, type a unit for the factor if applicable.

The Factor Unit is only a text label; however, it must start with a letter.

- 4. For each numeric value that you want to add to the Values list, do the following:
 - a. In the box next to the Add button, type a numeric value.

The Add button becomes available (Figure 39).

Figure 39.	Entering	numeric va	alues ir	the va	lues hox
riguie 55.	Linconing	numene ve			1003 001

Replicate	Apply Cancel ×
Factor Unit:	
Values:	
	1
2 Add	Delete

b. Click Add.

The value appears in the Values list in ascending order.

5. To save the study factor, click **Apply**.

The factor editor collapses to show only the study factor name and the Values list (Figure 40).

Figure 40. Numeric factor with a list of values

Replicate	Edit 🗙
	1
	2

Adding Biological Replicate Study Factors

You can add study factors when you create a study, when you add files to an existing study, and on the Study Definition page of an existing study.

You can add only one biological replicate factor to a study. Use the biological replicate factor for nested statistical models—that is, for studies that include one study factor nested within another study factor.

To add a biological replicate factor

 From the menu bar at the top right of the Study Factors pane, choose Add > Biological Replicate Factor.

The biological replicate factor editor appears with [new factor] automatically selected.

Figure 41. Biological replicate factor editor

8 [new factor]	Apply Cancel 🗙
Items:	
	Add Delete

2. Type a factor name to replace [new factor], for example, Rat.

- 3. For each item that you want to add to the Items list, do the following:
 - a. In the Items box (next to the Add button), type a study factor item, for example, **Rat 1**.

The Add button becomes available.

Figure 42. Biological replicate factor editor with added items

8 8 8	Rat	Apply Cancel 🗙
Items		
		Rat 1
		Rat 2
		Rat 3

b. Click Add.

The current item appears in the Items list.

4. To save the study factor, click **Apply**.

The factor editor collapses to show only the study factor name and the Items list. The items appear in alphabetical order.

Deleting Study Factors

To delete a study factor

1. In the Study Factors pane, click \mathbf{X} in the title bar of the factor.

Note Because the application cannot recognize whether a study factor is in use, this prompt appears even when you attempt to delete an undefined study factor.

2. At the prompt, click **Yes** to delete the study factor.

Duplicating Study Factors

To create a new study factor by using the Copy and Paste commands

1. In the Study Factors pane, select the factor that you want to copy.

The title bar of the selected factor turns blue.

- 2. Click Copy.
- 3. Click Paste.

A copy of the selected factor appears.

Editing Study Factors

You can edit study factors as you create a new study with the wizard, in Study Factors area of the Input File Characterization dialog box when you add more files to an existing study, and in the Study Factors area on the Study Definition page of an existing study.

To edit a study factor

1. In the factor title bar, click **Edit**.

The text entry box and the Add and Delete buttons appear. For a numeric factor, the Factor Unit box also appears.

- 2. To change the unit for a numeric factor, select the current unit and type a new unit.
- 3. To add more entries to the Items or Values list, type alphanumeric text in the appropriate box, and then click **Add**.
- 4. To delete an entry, select the entry and click **Delete**.

When an entry is in use, it is unavailable. To delete a value that is in use, you must first undo its assignment to any sample.

Automatically Assigning Study Factor Values

After you set up the study factors for a study, you can assign the study factor values to each sample. If the input file names follow a consistent pattern and the study factor values are completely defined, clicking Assign on the Input File Characterization page (or dialog box) assigns the study factor values to the samples.

Note You can only manually assign study factor values on the Samples page of an existing study.

- To automatically assign the study factor values to a sample set
- 1. If you have not already set up the study factors, set them up.
- 2. In the command bar of the Input File Characterization page or dialog box, click Assign.
- 3. Check the study factor columns, verify the sample assignments, and manually assign the study factor values if necessary.

Manually Selecting the Study Factor Values

After you set up the study factors for the study, you can select the study factor values for each sample. If the input file names do not follow a consistent pattern or do not include the study factor values or you are selecting the values on the Samples page of an existing study, you have to manually select these values.

✤ To manually select the study factor values

On the Input File Characterization page (or dialog box) or on the Samples page of an existing study, do the following:

- To select the factor values for a single sample, select the appropriate value from the list in each factor column.
- To select the same value for a consecutive sample range, drag the pointer across the rows of interest. Then right-click and choose **Set** *Factor* **To** > *Value*.
- To select the same value for nonconsecutive samples, hold down the CTRL key and click the samples of interest. Then right-click and choose **Set** *Factor* **To** > *Value*.

Resetting the Sample Assignments

On the Input File Characterization page (or dialog box), you can use the Reset button in the command bar to automatically clear the sample assignments. The Samples page of a study does not have a command bar.

* To clear the assignments in the Samples pane

In the command bar, click Reset.

Clicking Reset resets the Sample Type and study factor assignments—that is, it resets the sample type to Sample and the study factor values to n/a for all the samples.

Note To edit the values for a study factor, you must first clear the sample assignments if the values are assigned to samples.

Selecting the Sample Types

You can select the sample type for the study samples from any of these locations:

- The Samples pane of the wizard's Input File Characterization page
- The Samples pane of the Input File Characterization dialog box
- The Samples page of an existing study

By default, the selected sample type for each input file is Sample. Clicking Assign on the Input File Characterization page (or dialog box) automatically assigns the Blank sample type to samples with a file name that includes "blank" as a delimited value (for example, solvent_blank_1.raw, where the underscore character is the delimiter).

* To select a sample type other than the default assignment

In the Samples pane (any location), do one of the following:

• To select the sample type for a single sample, in the Sample Type column, select Sample, Control, Blank, Quality Control, Identification Only, Standard, or Labeled from the list.

San	nple	-
Sar	mple	
Co	ntrol	
Bla	nk	
Qu	ality Con	ntrol
Ide	ntificatio	on Only
Sta	ndard	
Lał	peled	

- To assign the same sample type to a consecutive sample range, use the SHIFT key to select a range of samples. Then, right-click and choose **Set Sample Type To** > *Sample Type*.
- To assign the same sample type to nonconsecutive samples, use the CTRL key to select the samples. Then, right-click and choose **Set Sample Type To** > *Sample Type*.

Tip To select a row, click a column without a list.

Related Topics

• Sample Types

Using the Regular Expression Builder to Extract the Study Factors

Use the Extract Sample Information From Sample Names dialog box to automate the setup and assignment of the study factors. This dialog box is accessible from the Input File Characterization page of the New Study and Analysis Wizard or from the Input File Characterization dialog box that opens when you add input files to a study.

Note For the extraction and assignment process to work, you must select the study factor portions of an example input file name and define these portions appropriately as categorical, numerical, or biological replicate factors.

When the undefined portions of the file name are not exactly the same for all the samples that you want to characterize, you can mark the text to be ignored or you can manually assign or edit the study factor values for these samples after returning to the Input File Characterization page or dialog box.

* To extract the sample information from the sample names

- 1. Open the Input File Characterization page (step 4 of 6) of the wizard or the Input File Characterization dialog box from within an existing study.
- 2. Click Advanced.

The Extract Sample Information From Sample Names dialog box opens with the first file name displayed in the Regular Expression Builder box and the file names of all the selected input files listed below it (Figure 43).



Extract Sample Information From Sample N Set up and assign the study factors, using the			
For information about regular expression op- click Help or press the F1 key.			Inserts a regular expression
Regular Expression Builder [0	-9] [A-Z] [/	A-Z] Text	that does the following:
Blank_01			 Ignores a text string. Defines a biological variable.
Sample Name			 Defines a categorical variable.
Blank_01			• Defines a numerical variable.
Blank_02			
Urine_0-4hr_01			
Urine_0-4hr_02			
Urine_4-8hr_01			
Urine_4-8hr_02			
Urine_8-12hr_01			
Urine_8-12hr_02			
Urine_Predose_01 Urine_Predose_02			
Unine_Predose_02			
<i>7</i>	ОК	Cancel	

3. If the Regular Expression Builder box does not contain a representative sample name, right-click a representative sample name in the Sample Name column and choose **Use** Name of Selected Sample as Regular Expression Template (Figure 44).

Figure 44. Urine_0-4hr_01 sample name selected

② Extract Sample Infor	rmation From Sample Names		×
	study factors, using the Regular E t regular expression operators and	xpression Builder. I examples of their use, click Help or press the F1 k	key.
Regular Expression B	uilder	[0-9] [A-Z] [A-Z] Text	
Blank_01			
Sample Name			
Blank_01			
Blank_02			
Urine_0-4hr_01			
Urine_0-4hr_02	Use Name of Selected Sa	ample as Regular Expression Template.	
Urine_4-8hr_01		45	-

The selected sample name replaces the nonrepresentative sample name (Figure 45).

Figure 45. Urine 0-4hr file name to be used as the regular expression template

Ŋ	Extract Sample Information From Sample Names
S	et up and assign the study factors, using the Regular Expression Builder.
F	or information about regular expression operators and examples of their use, click Help or press the F1 key.
1	Regular Expression Builder [0-9] [A-Z] [A-Z] Text
	Urine_0-4hr_01

- 4. For a representative sample name, do the following:
 - For each categorical study factor, select the corresponding study factor item and click [A-Z].

For example, select **0-4hr** and click **[A-Z]**.

The CategoricalVariable_1 column appears to the right of the sample names. In this example, the application extracts the categorical study factor for the

Urine_Predose_01 sample. The column is not populated with study factor values for the remaining samples for two reasons:

- The remaining portions of the sample names are not the same (for example, _01 and _02).
- The default expression for a categorical factor does not recognize the hyphen special character in the 0-4hr, 4-8hr, and 8-12hr time points.
- For each numeric study factor, select the corresponding study factor value and click **[0-9]**. For example, select **1** and click **[0-9]**.

For this example, the following expression replaces the file name:

```
Urine_(?<CategoricalVariable_1>[0-9A-Z]+)_ (?<NumericalVariable_1>[0-9]+)
```

-and-

Only the study factor values for these file names appear in the table: Urine_Predose_01 and Urine_Predose_02 (Figure 46).

Figure 46 shows the new regular expression in the regular expression builder box. This expression cannot interpret the time points that include hyphens (0-4hr, 4-8hr, 8-12hr). Only "Predose" matches the expression, which is looking for an alphanumeric string.

Figure 46. Building a regular expression with the default categorical and numeric operators

Regular Expression Bui	der		[0-9] [A-Z]	[A-Z] Text
Urine_(? <categoricalvariable_1>[0-9A-Z]+)_0(?<numericalvariable_1>[0-9]+)</numericalvariable_1></categoricalvariable_1>				9]+)
Sample Name	CategoricalVariable_1 <	Numeric	NumericalVariable_1	▼ INumeric
Blank_01				
Blank_02				
Urine_0-4hr_01				
Urine_0-4hr_02				
Urine_4-8hr_01				
Urine_4-8hr_02				
Urine_8-12hr_01				
Urine_8-12hr_02				
Urine_Predose_01	Predose		1	
Urine_Predose_02	Predose		2	

5. If the study factor columns do not populate as expected, modify the regular expression operators within the square brackets.

Operator	Quantifier	Matching pattern
[0-9]	none	A single integer
[0-9]	+	One or more integers
[A-Z]	none	One alphabetic character
[0-9A-Z]	none	One alphanumeric character
[0-9A-Z]	+	One or more alphanumeric characters
[0-9A-Z-]	+	One or more alphanumeric characters, one or more hyphens, or both (for example, 555-0000)
[0-9A-Z] [0-9A-Z]	+	One or more alphanumeric characters, a space, and one or more alphanumeric characters (for example, ABC 123)

Table 10. Examples of regular expression operators

For example, to populate the rows for the other file names with defined time periods, add a hyphen to the operators in the categorical factor's square brackets. Figure 47 shows the hyphen character added to the operator set for a categorical factor.

Figure 47. Hyphen building block added to the regular expression for a categorical factor

Original operator set = [0-9A-Z] — Modified operator set = [0-9A-Z-]

Extract Sample Information F	rom Sample Names					
et up and assign the study fac or information about regular e				ick Help or pre	ess the F1 k	ey.
Regular Expression Builder			[0-	·9] [[A-Z]	[A-Z]	Text
Urine_(? <categoricalva< th=""><th></th><th>_</th><th></th><th></th><th></th><th></th></categoricalva<>		_				
Sample Name	Categorica	Variał 🔻	Numeric	Numerical	Variab 🔻	Numeric
Blank_01						
Blank_02						
Urine_0-4hr_01	0-4hr			1		
Urine_0-4hr_02	0-4hr			2		
Urine_4-8hr_01	4-8hr			1		
Urine_4-8hr_02	4-8hr	4-8hr		2		
Urine_8-12hr_01	8-12hr	8-12hr		1		
Urine_8-12hr_02	8-12hr			2		
Urine_Predose_01	Predose			1		
Urine_Predose_02	Predose			2		

 If the sample names include extra text that differs from sample to sample and does not define a study variable, exclude this text from the regular expression by selecting it and clicking Text.

Note Figure 48 and Figure 49 show a different sample set than the previous figures.

For example, in the sample set shown in Figure 48, exclude all the characters to the right of the categorical variable by selecting _01_MDF and clicking _____.

Figure 48. Selecting the text that you want to exclude from the expression

🔰 Extract Sample Information From Sample Names 🛛 🗖	
Set up and assign the study factors, using the Regular Expression For information about regular expression operators and examples	
Regular Expression Builder [0-9] [A-Z]	Text
Urine_0-3_01_MDF	regular expression snippet to ignore text pattern.
Sample Name Urine_0-3_01_MDF Urine_0-3_GSH_01 Urine_0-3_GSH_0bs_01 Urine_3-5_GSH_PhII_01 Urine_3-5_GSH_PhII_01 Urine_5-7_01 Urine_7-9_01	The application ignores the extra text when extracting the study factor values from these sample names.
ØK Can	cel

The regular expression builder replaces the selected text with the following expression:

(?:.+)?

Figure 49 shows the result of these actions:

- Selecting the time period (0-3) as a categorical variable
- Adding a hyphen to the categorical variable expression
- Selecting "_01_MDF" as text to ignore

Figure 49. Regular expression that defines the categorical variable and the text to ignore

or information about regular expression	operators and examples of their use, cli	ick Help or press the F1 key.
Regular Expression Builder [0-9] [A-Z] Test		
Urine_(? <categoricalvariable_1>[0-9A-Z-]+)(?:.+)?</categoricalvariable_1>		
Sample Name	CategoricalVariable_1	▼
Urine_0-3_01_MDF	0-3	
Urine_0-3_GSH_01	0-3	
Urine_0-3_GSH_Obs_01	0-3	
Urine_0-3_GSH_PhII_01	0-3	
Urine_3-5_01	3-5	
Urine_3-5_GSH_PhII_01	3-5	
Urine_5-7_01	5-7	
Urine_7-9_01	7-9	

7. To enter the study factor names, type the study factor names in the column-heading boxes.

For example, replace CategoricalVariable_1 with **Time Period** and NumericalVariable_1 with **Replicate** (Figure 47 and Figure 50).

Figure 50. Study factors renamed

Extract Sample Information From S	Sample Names	
Set up and assign the study factors, u For information about regular expres	using the Regular Expression Builder. ssion operators and examples of their u	use, click Help or press the F1 key.
Regular Expression Builder		[0-9] [A-Z] [A-Z] Text
Urine_(? <categoricalvariab< td=""><td>ole_1>[0-9A-Z-]+)_0(?<nume< td=""><td>ericalVariable_1>[0-9]+)</td></nume<></td></categoricalvariab<>	ole_1>[0-9A-Z-]+)_0(? <nume< td=""><td>ericalVariable_1>[0-9]+)</td></nume<>	ericalVariable_1>[0-9]+)
Sample Name	Time Period Num	neric Replicate 🔻 🗹 Numeric

Edited study factor names

Tip You can edit the study factor names in two ways:

- In the column-heading boxes, replace the default column headings with the study factor names.
- On the Input File Characterization page, edit the study factors.

8. Click **OK** to return to the Input File Characterization page where you can modify the values in the study variable columns if necessary.

Tip The Regular Expression Builder does not assign sample types. To assign Blank as the sample type, click Assign on the Input File Characterization page after selecting the appropriate delimiters.

For example, select the **Underscore** check box on the Input File Characterization page and click **Assign**, for the following file name: Solvent_Blank_1.

 Table 11 describes the parameters in the Extract Sample Information From Sample Names

 dialog box.

Table 11. Extract Sample Information From Sample Names dialog box parameters (Sheet 1 of 2)

Parameter	Description
-----------	-------------

Regular Expression Builder

Initially displays the first file name in the Sample Name list. Use this file name (or a different file name in the list) to build a regular expression that extracts the study factor values for the file names in the Sample Name list.

Buttons	
[0-9]	Identifies a numeric factor.
	Clicking [0-9] replaces the selected text with the following expression: (? <numericalvariable_1>[0-9]+)</numericalvariable_1>
[A-Z]	Identifies a categorical factor.
	Clicking [A-Z] replaces the selected text with the following expression: (? <categoricalvariable_1>[0-9A-Z]+)</categoricalvariable_1>
[A-Z]	Identifies a biological replicate factor.
	Clicking [A-Z] replaces the selected text with the following expression: (? <replicatevariable_1>[0-9A-Z]+)</replicatevariable_1>
Text	Excludes text that does not define a study variable and does not follow a pattern.
	Clicking Text replaces the selected text with the following expression: (?:.+)?
Columns	
Sample Name	Displays the file names of the selected input files.
CategoricalVariable_1	Displays the extracted items for this factor. You can edit the column heading.

Parameter	Description
NumericalVariable_1	Displays the extracted values for this factor. You can edit the column heading.
ReplicateVariable_1	Displays the extracted values for this factor. You can edit the column heading.

Table 11. Extract Sample Information From Sample Names dialog box parameters (Sheet 2 of 2)

Setting Up the Sample Groups and Ratios for a New Analysis

Use the Sample Groups and Ratios page (step 5 of 6) of the New Study and Analysis Wizard to set up the sample groups that you want to compare and the group ratios that you want to include in the result file.

Note You can also use the Sample Groups and Ratios page within an existing study to set up the sample groups and ratios. However, when you start a new analysis in an existing study, you cannot set up the sample groups without first adding the input files to the Files for Analysis area of the Analysis pane.

IMPORTANT Do not create ratios with sample groups that have an n/a assignment for a study variable. To use a sample group in a ratio, you must replace the n/a assignment with a study factor value.

Follow these topics as needed:

- Opening the Wizard's Sample Groups and Ratios Page
- Selecting the Study Variables for Sample Grouping
- Fixing Assignment Errors for Unset Study Variables
- Changing the Hierarchy of the Study Variables
- Individually Setting Up the Group Ratios
- Using the Bulk Ratio Generation Feature to Set Up the Group Ratios

Opening the Wizard's Sample Groups and Ratios Page

- * To open the Sample Groups and Ratios page of the wizard
- 1. From the menu bar, choose File > New Study and Analysis and click Next.
- 2. Make the appropriate selections on the Study Name and Processing Workflow page and click **Next**.

The Input File Selection page opens.

3. Add all the input files for the study or add the input files for the current analysis and click **Next**.

If you selected a study template when you created the study, the Study Factors area of the Input File Characterization page contains one or more factors. Set up and assign the study factors and select the sample type for each sample.

4. Click Next.

The Sample Groups and Ratios page opens. The Study Variables area contains a File check box and a Sample Type check box. If you added study factors on the Input File Characterization page, the Study Variables area also contains an additional check box for each study factor.

Figure 51 shows a Sample Groups and Ratios page with generated sample groups and ratios.

Sample Group and Ratio Specification	Generated Sample Groups
Study Variables	n/a *
File	Blank n/a 1 F1: Solvent_Blank_1
Time Period	Blank n/a 2 F2: Solvent_Blank_2
Replicate	0-4hr
Sample Type	Sample 0-4hr 1 F3: Urine_0-4hr_1
	Sample 0-4hr 2 F4: Urine_0-4hr_2
- Manual Ratio Generation	4-8hr
Numerator:	Sample 4-8hr 1 F5: Urine_4-8hr_1
Add Ratio	Sample 4-8hr 2 F6: Urine_4-8hr_2
Denominator:	8-12hr
Bulk Ratio Generation	Sample 8-12hr 1 F7: Urine_8-12hr_1
Denominators to be used:	Sample 8-12hr 2 F8: Urine_8-12hr_2
Time Period : 0-4hr	Predose
Time Period : 4-8hr Time Period : 8-12hr	Sample Predose 1 F9: Urine_Predose_1
Time Period : 8-1211	Sample Predose 2 F10: Urine_Predose_2
	Generated Ratios 🐹 Clear Al
	× 0-4hr / Predose
	X 4-8hr / Predose
Add Ratios	X 8-12hr / Predose

Figure 51. Sample Groups and Ratios page of the wizard

Selecting the Study Variables for Sample Grouping

You select the study variables that you want to group on the Grouping & Ratios page of an existing analysis or the Sample Groups and Ratios page of the New Study and Analysis wizard.

* To select the study variables for sample grouping

In the Study Variables area of the Sample Groups and Ratios page, select the study variable or variables for sample grouping.

Selected study variables have a light green background, a dark green handle ([]) on the left, and a Sorting icon ([]) on the right.

The Generated Sample Groups area displays the generated sample groups. The hierarchy of the study variables affects the sample group names and the denominator list for bulk ratio generation (see To change the hierarchy of the study variables).

The naming scheme is as follows:

- Group names (green) consist of the common values for the selected study variables.
- Sample names (blue) consist of a unique ID and the input file name.

Following data processing, the result file displays the chromatographic peak areas for individual samples and the named sample groups.

Fixing Assignment Errors for Unset Study Variables

You select the study variables that you want to group on the Grouping & Ratios page of an existing analysis or the Sample Groups and Ratios page of the New Study and Analysis wizard.

When selecting a study variable generates an N/A group with Sample, Control, or Standard sample types, the application highlights the selected study variable and the N/A group in red and displays an error message.

To fix an assignment error

- 1. Do one of the following:
 - In the New Study and Analysis Wizard, return to the Input File Characterization page.
 - From an existing study, open the Samples page.
- 2. Do one of the following:
 - Assign study factor values to the Sample, Control, or Standard sample types. If necessary, create new study factor values for these samples.
 - Change the sample type assignment for samples without a study factor value to **Blank** or **Identification Only**.

- 3. Do one of the following:
 - In the New Study and Analysis Wizard, return to the Sample Groups and Ratios page and verify the assignments.
 - In an existing study where you have started a new analysis, return to the Grouping & Ratios page.

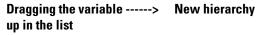
Changing the Hierarchy of the Study Variables

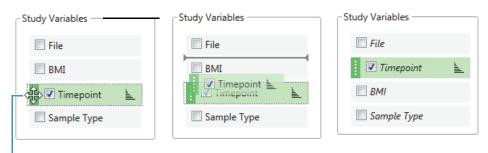
You can change the hierarchy of the study variables on the Grouping & Ratios page of an existing analysis or the Sample Groups and Ratios page of the New Study and Analysis wizard.

* To change the hierarchy of the study variables

Place the pointer over the handle (1) to the left of a variable name. When the move cursor (4) appears, drag the variable up or down in the list.

Original hierarchy ----->





Move cursor

Figure 52 and Figure 53 show the effect of changing the study variable hierarchy for a study where either the extraction solvent (IPA or WFI) or the o-ring type (E3609, FF350, or HF355) is the primary study variable.

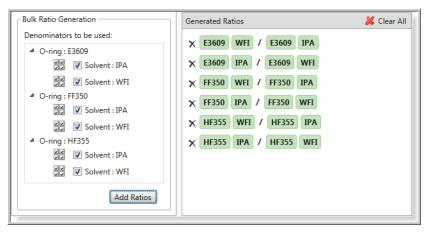
The ratios in Figure 52 (where extraction solvent is the primary variable) compare the compounds extracted with the same solvent from different o-ring types.

The ratios in Figure 53 (where o-ring type is the primary variable) compare the compounds extracted from each o-ring type with the two solvents.

Bulk Ratio Generation	Generated Ratios	样 Clear All
Denominators to be used:	× IPA FF350 / IPA E3609	
Solvent : IPA		
🖉 🗹 O-ring : E3609	X IPA HF355 / IPA E3609	
🖉 🔽 O-ring : FF350	X IPA E3609 / IPA FF350	
VV O-ring : HF355	× IPA HF355 / IPA FF350	
▲ Solvent : WFI		
🖉 🗹 O-ring : E3609	X IPA E3609 / IPA HF355	
🖉 🗹 O-ring : FF350	× IPA FF350 / IPA HF355	
V O-ring : HF355	X WFI FF350 / WFI E3609	
Add Ratios	X WFI HF355 / WFI E3609	
	X WFI E3609 / WFI FF350	
	X WFI HF355 / WFI FF350	
	X WFI E3609 / WFI HF355	
	X WFI FF350 / WFI HF355	

Figure 52. Extraction solvent is the primary variable

Figure 53. O-ring type is the primary study variable



Individually Setting Up the Group Ratios

You can set up the group ratios on the Grouping & Ratios page of an existing analysis or the Sample Groups and Ratios page of the New Study and Analysis wizard.

To set up group ratios one-by-one

In the Manual Ratio Generation area, select a group from the Numerator list, select a group from the Denominator list, and click **Add Ratio**. Repeat this step for each pair of groups that you want to compare.

Using the Bulk Ratio Generation Feature to Set Up the Group Ratios

You can set up the group ratios on the Grouping & Ratios page of an existing analysis or the Sample Groups and Ratios page of the New Study and Analysis wizard.

The sort order of the study variables affects the denominators to be used in the Bulk Ratio Generation area.

To automatically set up multiple ratios

1. To change the sort order of the study variables, click the **Sorting** icon for the study variable that you want to sort by and choose **Sort Ascending** or **Sort Descending**. If you do not want to apply a sort order, choose **No Sorting**.

Time Period	E	
		No Sorting
	1	Sort Ascending
	llin.	Sort Descending

2. In the Bulk Ratio Generation area, select one or more denominators and click Add Ratios.

When you select more than one variable in the Study Variables area and the selected variables have more than one value, the Select/Deselect Item in All Groups icon, the appears when you place the pointer near a denominator in the Denominators to Be Used area.

To automatically select multiple check boxes for the same value in all groups, select one of the check boxes for the value, and then click the **Select/Deselect Item in All Groups** icon,

- 3. To modify the group ratio list, do either of the following:
 - To remove a ratio, click the delete icon (X) to its left.
 - To clear the entire list, click **Clear All**.

Preparing to Submit a Run to the Job Queue

The first time you use the New Study and Analysis Wizard, read the instructions on the final page of the wizard.

Clicking Finish on the Input File Selection page (after selecting input files), on the Input File Characterization page, or on the final wizard page opens the study page and the Analysis pane. The Analysis pane contains the selected input files. If you selected a processing workflow on the Study Name and Processing page of the wizard, the Workflow box on the Analysis pane displays the name of the processing workflow, and the Workflows page contains the processing workflow.

Before starting a run, you can edit the study, the processing workflow, the input file list, and the result file name. Some of the defined processing workflows require customization. For example, for a targeted analysis, you must select the target compounds for the Generate Expected Compounds node.

If the processing workflow is valid and the Analysis pane contains one or more input files, the Run command at the top right of the Analysis pane is available.

If the Run command is unavailable, you must fix the analysis errors.

Related Topics

• Troubleshooting Common Analysis Errors

About Accessing the Wizard Help

The New Study and Analysis Wizard has two types of Help:

- Embedded instructions that appear to the left of each wizard page
- Context-sensitive Help topics that are part of the Help system

Task	Action
Display the embedded instructions for the current page.	In the lower left corner of the wizard, click the light bulb icon (\bigcirc).
Hide the embedded instructions for the current page.	In the lower left corner of the wizard, click the light bulb icon (🛐).
Open the context-sensitive Help topic for the current page.	Press the F1 key (F1) on your computer keyboard.

Editing Existing Studies

In the Compound Discoverer application, you process your raw data files (run analyses) within the study environment. These topics describe the study features and how to edit an existing Compound Discoverer study.

Contents

- About Existing Study Files
- Opening an Existing Study
- Study Page Commands and Tabs
- Editing the Study Factors and the Study Description
- Adding Input Files to an Existing Study
- Removing Input Files or Updating Their Location
- Editing the Sample Type and Study Factor Values
- Saving a Study File

About Existing Study Files

When you create a study with the New Study and Analysis Wizard, the application creates a study file (.cdStudy) and a study folder. The study file contains a list of input files (Xcalibur RAW files) with their associated sample information and a list of analyses with their associated result files (.cdResult). The sample information includes the sample type of each input file and the relationship between the input files.

When you open an existing study, it opens as a tabbed document with a command bar and a set of tabbed pages: Study Definition, Input Files, Samples, and Analysis Results.

The Input Files page of a study tracks the status of the input files. These input files typically reside outside the study folder, for example, on a shared server or on the data system computer where they were acquired. If you delete or rename an input file from the specified folder location after you add it to a study, this warning symbol, **()**, appears to the left of the ID column for the deleted file. Placing the cursor over the warning symbol displays the missing file's expected location and instructions for resolving the issue.

4

The Analysis Results page of a study tracks the result files generated by analyses run within the study. Result files reside within the study's folder.

Related Topics

• Resolving the Location of a Study's Input Files

Opening an Existing Study

You can open an existing study from the Start Page or the Compound Discoverer window.

To open an existing study

Do one of the following:

- From the Start Page, under What Would You Like to Do?, click **Open Study** to open the Open Study dialog box. Select a study file and click **Open**.
- From the Start Page, under Recent Studies, click the study name of interest.

The Start Page lists the 20 most recent studies. The study name appears in blue hypertext. Placing the cursor over a study name underlines it (Figure 54).



		Me		Study name n blue hypert	ext	Toolbar
Compound	d Discoverer 3.1.0					
<u>F</u> ile <u>R</u> eporti	ng <u>L</u> ists & Libraries <u>V</u> iew <u>W</u> indow	<u>H</u> elp				
	: <mark>■</mark>	E [] 8	: <u>(</u> 14 : °,	þ. °. : © C	*: 🛉 🖒 🖏	
	Compound vould you like to do?		SCOVEI Studies		1 Results	-
r <mark>i i</mark> n	New Study and Analysis Create a new study and analysis from scratch.	8 0	ZDF Study C:\Studies\ZDF Study	ø	Fatty vs Lean C:\Studies\ZDF Study	
æ	Open Study Open an existing study.		Omeprazole Stud C:\Studies\Omeprazole		Target Omeprazole C:\Studies\Omeprazole Stud	y
0	Open Result Open an existing result file.					

From the application menu bar, choose File > Open Study to open the Open Study dialog box. Then, select a study file and click Open.

-or-

• From the application toolbar, click the **Open an Existing Study** icon, ^{III}, to open the Open Study dialog box. Select a study file and click **Open**.

The study opens to the Analysis Results page. The study tab includes an image of two racked test tubes, f = 1, on the left, the study name in the middle, and a close icon, \mathbf{X} , on the right. When you make changes to a study, an asterisk (*) appears to the right of the study name to indicate unsaved changes (Figure 55).

Tip See the option for "Turning Off the Auto-Save Feature for Studies" on page 567.

Figure 55. Analysis Results page of a study

1	Start Pa	age × III ZDF Study* ×					
🙀 Add Files 💥 Remove Files 🔍 Open Containing Folder New Analysis 🍯 Open Analysis Template							
Stud	y Defin	ition Input Files Samples	Analysis Results				
6	Open R	esult 🛯 🎲 Reprocess					
Error	ID 🔺	Execution State	Creation Date	File Name	File Type	File Size	Description
		•	II	I .			
	1	Completed (with Warnings)	5/18/2018 12:47 PM	Fatty vs Lean.cdResult	.cdResult	995,208 KB	Untargeted - Performs

Study Page Commands and Tabs

The study page includes a command bar and four tabbed pages (Table 12).

 Table 12.
 Study page parameters (Sheet 1 of 2)

Command or page	Description	
Commands		

The Add Files, New Analysis, and Open Analysis Template commands are independent of the active page within the study. The Remove Files command is only active when a file is selected on the Input Files page or the Analysis Results page.

Add Files	Opens the Open dialog box where you select the input files (Xcalibur RAW files) that you want to include in the study.
Remove Files	Executes one of two actions:
	• Removes the selected files on the Input Files page when it is the active page.
	• Removes the selected analysis results on the Analysis Results page when it is the active page.
Open Containing Folder	Opens the folder that contains the selected file.
	This command is only available for the Input Files and Analysis Results pages.

tab set on the study page. The Workflows page itself is emp This command is unavailable when the Analysis pane is opOpen Analysis TemplateOpens the Analysis Template dialog box where you select a analysis template. This command is unavailable when the Analysis pane is op Tabbed pages Use these pages as follows.Study DefinitionSet up study factors and view the study name, file location creation date, modification date, and description on this page lists the file ID, file name, file type, and sample information for each input file. Clicking Add Files opens the Input File Characterization dialog box and the Input Files page, if it is not already opeSamplesSelect the sample type and study factors for each raw data Access the result files created within the study, review the	Command or page	Description
Open Analysis TemplateOpens the Analysis Template dialog box where you select a analysis template. This command is unavailable when the Analysis pane is opTabbed pagesUse these pages as follows.Study DefinitionSet up study factors and view the study name, file location creation date, modification date, and description on this pa Input FilesInput FilesTrack the status and resolve the location of input files, as the page lists the file ID, file name, file type, and sample information for each input file.Clicking Add Files opens the Input File Characterization dialog box and the Input Files page, if it is not already opeSamplesAnalysis ResultsAccess the result files created within the study, review the	New Analysis	Opens the Analysis pane and adds the Workflows tab to the tab set on the study page. The Workflows page itself is empty.
Image: The second sec		This command is unavailable when the Analysis pane is open.
Tabbed pagesUse these pages as follows.Study DefinitionSet up study factors and view the study name, file location creation date, modification date, and description on this pageInput FilesTrack the status and resolve the location of input files, as the page lists the file ID, file name, file type, and sample information for each input file.Clicking Add Files opens the Input File Characterization dialog box and the Input Files page, if it is not already opeSamplesSelect the sample type and study factors for each raw data Access the result files created within the study, review the	Open Analysis Template	Opens the Analysis Template dialog box where you select an analysis template.
Use these pages as follows. Study Definition Set up study factors and view the study name, file location creation date, modification date, and description on this pa- Input Files Track the status and resolve the location of input files, as the page lists the file ID, file name, file type, and sample information for each input file. Clicking Add Files opens the Input File Characterization dialog box and the Input Files page, if it is not already ope Samples Select the sample type and study factors for each raw data and Analysis Results Access the result files created within the study, review the		This command is unavailable when the Analysis pane is open
Study DefinitionSet up study factors and view the study name, file location creation date, modification date, and description on this paInput FilesTrack the status and resolve the location of input files, as the page lists the file ID, file name, file type, and sample information for each input file.Clicking Add Files opens the Input File Characterization dialog box and the Input Files page, if it is not already opeSamplesSelect the sample type and study factors for each raw data Analysis Results	Tabbed pages	
creation date, modification date, and description on this particularInput FilesTrack the status and resolve the location of input files, as the page lists the file ID, file name, file type, and sample information for each input file.Clicking Add Files opens the Input File Characterization dialog box and the Input Files page, if it is not already opeSamplesSelect the sample type and study factors for each raw dataAnalysis ResultsAccess the result files created within the study, review the	Use these pages as follows.	
Page lists the file ID, file name, file type, and sample information for each input file.Clicking Add Files opens the Input File Characterization dialog box and the Input Files page, if it is not already opeSamplesSelect the sample type and study factors for each raw data Analysis ResultsAccess the result files created within the study, review the	Study Definition	Set up study factors and view the study name, file location, creation date, modification date, and description on this page
dialog box and the Input Files page, if it is not already opeSamplesSelect the sample type and study factors for each raw dataAnalysis ResultsAccess the result files created within the study, review the	Input Files	
Analysis Results Access the result files created within the study, review the		Clicking Add Files opens the Input File Characterization dialog box and the Input Files page, if it is not already open.
	Samples	Select the sample type and study factors for each raw data file
analysis details for the results mes, and reprocess an analysis	Analysis Results	Access the result files created within the study, review the analysis details for the results files, and reprocess an analysis.

Table 12. Study page parameters (Sheet 2 of 2)

Related Topics

- Removing Input Files or Updating Their Location
- Editing the Study Factors and the Study Description
- Editing the Sample Type and Study Factor Values
- Reviewing or Reprocessing an Analysis

Editing the Study Factors and the Study Description

Use the Study Definition page of a study to edit the existing study factors, to set up new study factors, and to edit or add a description of the study.

Follow these topics as needed:

- Editing the Study Description
- Adding, Deleting, or Editing Study Factors in an Existing Study
- Study Definition Page of a Study

Note If you selected a study template with study factors when you created the study or added study factors by using the Input File Characterization page of the New Study and Analysis wizard, the Study Factors area contains the defined factors.

Editing the Study Description

You can edit the description of a study on its Study Definition page.

✤ To add or edit a description of the study

Type or edit the description in the Study description area.

Adding, Deleting, or Editing Study Factors in an Existing Study

To add, delete, or edit study factors in an existing study, see the following subtopics in "Characterizing New Input Files" on page 77.

- Adding Numeric Study Factors
- Adding Biological Replicate Study Factors
- Deleting Study Factors
- Editing Study Factors
- Duplicating Study Factors

Study Definition Page of a Study

Table 13 describes the parameters on the Study Definition page.

Table 13. Study Definition page parameters (Sheet 1 of 3)

Parameter	Description
Study Summary pane	
Study Name	Displays the study name.

Parameter	Description
Study Directory	Displays the file location where the study is stored.
Last Changed	Displays the date and time of the last saved change to the study.
Creation Date	Displays the creation date of the study file.
Study Description pane	
Use this pane to enter and st	ore a description of the current study.
Study Factors pane	
Use this pane to set up or ed	it the study factors.
Menu commands	
Paste	Pastes the entries in the copied factors below the existing factors.
Сору	Copies the selected factors to the Clipboard.
Add > Biological Replicate Factor	Opens a blank biological replicate editor. You can use the biological replicate factor to create nested studies.
Add > Categorical Factor	Opens a blank categorical factor editor.
Add > Numeric Factor	Opens a blank numeric factor editor. The numeric factor editor only accepts numeric values.
Factor box	
Title bar	Displays the editable factor name.
Buttons and icons	
Apply	Saves the current entries in the factor editor.
Cancel	Closes the item or value entry box and removes any entries made during the current editing session. Does not remove previously saved entries.
Х	Deletes the factor from the study.
Add	Adds an item to a categorical factor or a numeric value to a numeric factor.
Delete	Deletes the selected item or value from the list in the respective Items or Values area.
Text entry boxes	
[new factor]	Type a factor name in this box.

Table 13. Study Definition page parameters (Sheet 2 of 3)

Parameter	Description
Item box	Type the name of an item that you want to add to the Items list for a categorical factor in this box.
Value box	Type a numeric value that you want to add to the Values list for a numeric factor in this box.

Table 13. Study Definition page parameters (Sheet 3 of 3)

Adding Input Files to an Existing Study

Use the Add Files button on the study command bar to add files to a study.

✤ To add input files to an existing study

- 1. Open the study of interest.
- 2. On the study command bar, click Add Files.
- 3. In the Add Files dialog box, select the files of interest and click Open.

The Input File Characterization dialog box opens.

- 4. Select the sample types and study factor values for the new samples.
- 5. Click OK.

The Input Files page of the study opens.

Related Topics

• Characterizing New Input Files

Removing Input Files or Updating Their Location

Use the Input Files page to remove or update the location of the input files in an existing study. The input files for the Compound Discoverer application are Xcalibur RAW files acquired by a Thermo Scientific HRAM mass spectrometer.

Tip To add input files to an analysis, drag the files of interest from the Input Files page to the Analysis pane (see "Setting Up a New Analysis from Within an Existing Study" on page 115).

Follow these topics as needed:

- Displaying the Location of a Study's Input Files
- Removing Input Files from a Study
- Resolving the Location of a Study's Input Files
- Input Files Page Parameters

Displaying the Location of a Study's Input Files

To display the file path for an input file

- 1. On a study's Input Files page, select the input file of interest.
- 2. Below the table, click **Show Details** (Figure 56).

Figure 56. Input Files page

Stud	ly Defin	ition Input Files Samples An	alysis Results				
rror	ID 🔺	Name	File Type	Sample Information			
•		•		II. •			
	F1	blank	.raw	Sample Type: [Blank], Phenotype: [n/a]			
	F2	Pooled_ddMS2	.raw	Sample Type: [Identification Only], Phenotype: [n/a]			
	F3	ZDF_Fatty_1	.raw	Sample Type: [Sample], Phenotype: [Fatty]			
	F4	ZDF_Fatty_2	.raw	Sample Type: [Sample], Phenotype: [Fatty]			
	F5	ZDF_Fatty_3	.raw	Sample Type: [Sample], Phenotype: [Fatty]			
	F6	ZDF_Lean_1	.raw	Sample Type: [Sample], Phenotype: [Lean]			
	F7	ZDF_Lean_2	.raw	Sample Type: [Sample], Phenotype: [Lean]			
	F8 ZDF_Lean_3 .raw Sample Type: [Sample], Phenotype: [Lean]						

Show/Hide Details command

The Samples page opens below the input files table.

3. Click the **Files** tab.

Figure 57 shows the hidden Files page that lists the selected input file's full file name, including its directory location.

Figure 57. File path of selected input file

	e Details es Files				
	File Name	File Size [kB	File Time		
	II. •			*	
F1.1	C:\Metabolomics Study\ZDF Study\blank.raw	347979	9/29/2015 2:3	8 PM	

Removing Input Files from a Study

To remove input files from a study

1. On the Input Files page, select the rows to remove and click **Remove Files** in the study command bar.

Depending on whether you have run an analysis with the selected input files, one of the following confirmation boxes appears:

- If you have not run an analysis, the Remove Input File confirmation box appears.
- If you have run an analysis, the Remove Analysis Result Files confirmation box appears.
- 2. Do one of the following:
 - In the Remove Input File confirmation box, click **Yes** to remove the input files from the study.

-or-

• In the Remove Analysis Result Files message box, click **Remove Files** to remove the input files and their associated analyses from the study.

When you remove an input file from the study, the analyses associated with the input file disappear from the Analysis Results page, but the result files (.cdResult) remain in the study folder.

Resolving the Location of a Study's Input Files

• To resolve the input files list when you move a study or the Xcalibur RAW files

Note If the network or directory path changes between the study and the raw data files, an exclamation symbol (**1**) appears to the left of the file ID.

- 1. On the study's Input Files page, display the details for the missing files and check their original location.
- 2. If you know where the files are currently stored, add the files to the study by using the Add Files command in the Study page command bar or by using the Find Missing Files command from the shortcut menu.

The Adding Files confirmation box opens. The progress remains at 0.0% until you click OK (Figure 58).

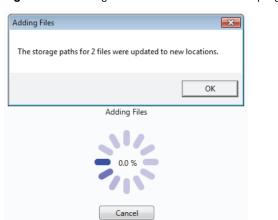


Figure 58. Adding Files confirmation box with progress information

3. Click **OK** to continue.

The Adding Duplicate Files message box opens. The progress again remains at 0.0% until you click OK.

4. Click **OK** to continue.

The application restores the connection between the study and the raw data files.

Input Files Page Parameters

Table 14 describes the parameters on the Input Files page.

Table 14. Input Files page parameters (Sheet 1 of 2)

Parameter or feature	Description
Show Details/Hide Details	Displays or hides the Samples and Files subpages.
Columns	
Error	Displays an exclamation mark, () , if the file is missing.
ID	Displays a unique ID in the following format: F#, where # is a unique integer. If you remove a file, and then add it again, the application updates the ID number.
Name	Displays the file name of the raw data file.
File Type	Displays the file type of the input file. The Compound Discoverer application supports Xcalibur RAW files (.raw).
Sample Information	Displays the sample type and any other selected study factors.
Hidden pages (Samples and Fil	es)
Samples page	
Sample	Displays a unique identifier for the input file.

Parameter or feature	Description
Sample Identifier	Displays the file name of the raw data file.
Sample Type	Displays the sample type (see "Sample Types" on page 68).
Study factor columns	Displays the study factor values. You can modify the study factor selections.
	Changing the factor values and sample types on the Input Files page also updates these items on the main Samples page.
Files page	
ID	Displays a unique ID (reserved for future implementation).
Name	Displays the file name and directory location of the raw data file.
Date Modified	Displays the acquisition date and time of the raw data file.
Size	Displays the size of the raw data file in bytes.
Shortcut menu for the Input Fi	les page
Set As Input File	Adds the selected input file to the Files for Analysis area in the Analysis pane.
	Available when the Analysis pane is open.
Find Missing Files	Opens the Browse for Folder dialog box. Browse to the folder where the file is stored and click OK . Then, click OK to close the confirmation message.
	Available when the Input Files table contains missing files.

Table 14. Input Files page parameters (Sheet 2 of 2)

Editing the Sample Type and Study Factor Values

If you have not already characterized the study samples or you want to change their characteristics, use the Samples page (Figure 59) of an existing study. To open a study's Samples page, click the Samples tab.

Figure 59. Samples page and its shortcut menu

	Stu	dy Definiti	on I	nput Files Samples Analysis Results				
	Erro	Samp 🔺	File	Sample Identifier	Sample Type	Ph	henotype	
ter				•	• •		J +	Copy With Headers Ctrl+C
N		S1	F1	blank	Blank •	Le	.ean 🔹	Сору
		S2	F2	Pooled_ddMS2	Identificatior *	n	n/a ▪	
		S3	F3	ZDF_Fatty_1	Sample *	Fa	atty 🔹	Clear Selection
		S4	F4	ZDF_Fatty_2	Sample +	Fa	atty 🔹	Cell Selection Mode
		S5	F5	ZDF_Fatty_3	Sample +	Fa	atty 🔹	Enable Row Grouping
		S6	F6	ZDF_Lean_1	Sample +	Le	.ean 🔹	
		S7	F7	ZDF_Lean_2	Sample +	Le	.ean 🔹	Set Sample Type to
		S8	F8	ZDF_Lean_3	Sample +	L	.ean 🔹	Set Phenotype to
	•	Show Asso	ciated	File				Set as Input File

To select or modify the sample types and study factor values or view the file information on the Samples page, follow these topics as needed:

- Editing the Sample Identifier on the Samples Page
- Editing the Sample Types on the Samples Page
- Editing the Study Factor Values on the Samples Page
- Viewing the File Information for Study Samples on the Samples Page

Editing the Sample Identifier on the Samples Page

To make it easier to identify each sample, you can modify the text in the Sample Identifier column.

* To edit the text in the Sample Identifier column

1. Point to the sample cell.

The Edit icon appears.

- 2. Click the **Edit** icon, **2**.
- 3. Place the cursor in the cell, and then type or paste the new text.

Editing the Sample Types on the Samples Page

* To edit the sample types

Tip To select a row, click a column that does not include a dropdown list.

Do any of the following:

- To select the sample type for a single sample, in the Sample Type column, select Sample, Control, Blank, Quality Control, Identification Only, Standard, or Labeled from the list.
- To assign the same sample type to a consecutive sample range, use the SHIFT key to select a range of samples. Then, right-click and choose **Set Sample Type To** > *Sample Type*.
- To assign the same sample type to nonconsecutive samples, use the CTRL key to select the samples. Then, right-click and choose **Set Sample Type**.

Related Topics

• Sample Types

Editing the Study Factor Values on the Samples Page

To edit the study factor values

Do any of the following:

- To select the factor values for a single sample, select the appropriate value from the list in each factor column.
- To select the same value for a consecutive sample range, drag the pointer across the rows of interest. Then right-click and choose **Set** *Factor* **To** > *Value*.
- To select the same value for nonconsecutive samples, hold down the CTRL key and click the samples of interest. Then right-click and choose **Set** *Factor* **To** > *Value*.

Tip To select a row, click a column that does not include a dropdown list.

Viewing the File Information for Study Samples on the Samples Page

* To view the file name and location for a selected sample

Click Show Associated File at the bottom left of the Samples page.

The following details appear—the full file name and location, the file size in kilobytes, and the acquisition time of the Xcalibur RAW file.

Samples Page Parameters

Table 15 describes the parameters on the Samples page.

 Table 15.
 Samples page parameters

Description
Shows or hides the file information for a selected sample.
Displays an error symbol.
Displays a unique number for the sample (S#).
Displays a unique number for the input file (F#).
Displays the file name of the raw data file.
Specifies the sample type for each sample.
Selections: Sample, Control, Blank, Quality Control, Identification Only, Standard, or Labeled.
Specifies the study factor values for each sample.
You can edit the study factor values.
Displays a unique file identifier.
Displays the file name and directory location of the Xcalibur RAW data file.
Displays the size of the Xcalibur RAW data file in bytes.
Displays the acquisition date and time of the Xcalibur RAW data file.

Snortcut menu commands

For information about the Copy with Headers, Copy, Clear Selection, and Cell Selection Mode commands, see "Working with the Application Tables" on page 579.

Set Sample Type To	Assigns the selected sample type to a sample range.
Set Factor To	Assigns the selected study factor value to a sample range.
Set As Input File	Adds the selected input file to the Files for Analysis area in the Analysis pane.
	Available when the Analysis pane is open.

Saving a Study File

By default, the application saves the study file when you close it. However, you can change this behavior by turning off the auto save feature.

IMPORTANT If you turn off the auto save feature, the application prompts you to save your changes when you attempt to close a study file with unsaved changes.

Unsaved changes include, for example, the last completed run on the Analysis Results page, new input files on the Input Files page, new study factors or study factor values on the Study Definition page, new sample assignments on the Samples page, and so on.

✤ To save a study file

Do one of the following:

From the menu bar, choose File > Save All, or in the toolbar, click the Save All Open Items icon,

-or-

Click the study tab to make it the active page. Then, from the menu bar, choose File
 > Save, or in the toolbar, click the Save the Currently Active Item icon,

Related Topics

• Turning Off the Auto-Save Feature for Studies

4 Editing Existing Studies Saving a Study File

5

Creating, Running, and Reprocessing Analyses

To set up and submit new analyses, review the details of completed analyses, or reprocess analyses, see these topics.

Contents

- Setting Up a New Analysis from Within an Existing Study
- Troubleshooting Common Analysis Errors
- Common Analysis Errors
- Submitting an Analysis to the Job Queue
- Common Validation Issues
- Working with the Job Queue
- Reviewing or Reprocessing an Analysis
- Analysis Pane Parameters

Setting Up a New Analysis from Within an Existing Study

You can set up a new analysis by beginning with an empty Workflows page, an empty Grouping & Ratios page, and an empty Analysis pane. Or, you can set up a new analysis by reprocessing an analysis from the Analysis Results page.

Tip You cannot set up the sample groups and ratios on the Grouping and Ratios page without first adding the appropriate input files to the Analysis pane.

To set up a new analysis from within an existing study, follow these topics in order:

- 1. Starting a New Analysis from Within an Existing Study
- 2. Selecting the Input Files for a New Analysis
- 3. Naming the Result File for a New Analysis
- 4. Setting Up the Sample Groups and Ratios for a New Analysis

For information about reprocessing an analysis, see "Reprocessing an Analysis" on page 131.

Starting a New Analysis from Within an Existing Study

This topic describes how to start a new analysis from within an existing study. For information about reprocessing an existing analysis, see "Reprocessing an Analysis" on page 131.

* To start a new analysis from within an existing study

- 1. Open the study.
- 2. In the study command bar, click New Analysis.

The Workflows tab and the Grouping & Ratios tab appear to the right of the study page tabs and an empty Analysis pane opens to the right of the tabbed pages. Both the Workflows and Grouping & Ratios pages are empty.

Note You can also start a new analysis by opening an analysis template.

- 3. To select or create a processing workflow for the analysis, do the following:
 - a. Click the Workflows tab.

<u></u>	ouping & Rat			I Save Common	Muto Layout 🛛 💥 Clea		Command bar
	rkflow:	open common	IN COVE	ing save common		_	
Des	cription:					* *	Workflow box
Wor	rkflow Tree						

- b. In the Workflow Tree area, create a new processing workflow or in the Workflows command bar, do one of the following to select an existing processing workflow:
 - To select a standard processing workflow from the Common Templates folder, click **Open Common**, click **Workflow Templates**, click the folder for your field of interest, select one of the provided processing workflow templates, and click **Open**.

The file name appears in the Workflow box, and a description of the processing workflow appears in the Description box.

• To select a processing workflow from another folder, click **Open**, locate the processing workflow file (.cdProcessingWF) of interest, and click **Open**.

The file name appears in the Workflow box.

-or-

• To load the processing workflow that you previously used to process a specific result file, click **Open**, locate the result file (.cdResult) of interest, and click **Open**.

4. Customize the processing workflow as needed.

Table 16 lists the workflow nodes that require custom settings.

Table 16. Workflow nodes that require customization

Node	Required customization					
Generate Expected Compounds	From the Compound list, select library compounds.					
Compounds	In the defined workflow templates, this selection is empty.					
Create FISh Trace	From the Compound list, select a library compound.					
	In the defined workflow templates, this selection is empty.					
Create Pattern Trace	In the Isotope Ratios box, define the ratio.					
	In the defined workflow templates, the ratio has been set to C15S.					
Search Mass Lists	From the Input Files list, select mass lists.					
	In the defined workflow templates, one of the preinstalled mass lists has been selected, according to the vertical market.					
Search mzVault	From the mzVault Library list, select mzVault libraries.					
	In the defined workflow templates, the preinstalled mzVault May 2019 library has been selected.					
Compound Class Scoring	From the Compound Classes list, select a Compound Class file.					
Pattern Scoring	For the Isotope Patterns parameter, set up an isotope pattern.					
	In the defined workflow templates, the pattern has been set to C15S. You can define and add more ratios.					

Thermo Fisher Scientific has optimized most of the parameter settings in the defined processing workflow templates by the area of study (vertical market). Table 17 lists the parameters that usually require a different setting, regardless of whether you use one of the defined templates or create your own processing workflow.

 Table 17. Parameter modifications for workflow nodes (Sheet 1 of 2)

Node	Parameter settings to optimize or modify
Create Analog Trace	Select the trace of interest.
Create Mass Trace	Select the trace of interest.
Generate Expected Compounds	Set up the dealkylation and transformation steps and select the appropriate ions.

Node	Parameter settings to optimize or modify
Search nodes	Select the libraries or lists of interest.
Detect Compounds	For the recommended minimum peak intensity settings for a specific mass spectrometer model, see Table 18.

Table 17. Parameter modifications for workflow nodes (Sheet 2 of 2)	Table 17.	Parameter	modifications	for workflow	nodes (Sheet 2 of 2)
---	-----------	-----------	---------------	--------------	----------------------

Table 18 lists the recommended range for the minimum peak intensity parameter. The optimal setting depends on the sensitivity of the mass spectrometer.

Table 18. Recommended minimum peak intensity range

Mass spectrometer	Minimum peak intensity (chromatographic peak height)
Q Exactive, Q Exactive Plus™, Q Exactive HF	500 000 to 1 000 000
Orbitrap Fusion, Orbitrap Lumos™, Orbitrap ID-X™	50 000 to 100 000
Exactive, Exactive Plus™, Orbitrap Elite™, Orbitrap Velos Pro™	100 000 to 500 000
LTQ Orbitrap XL™, LTQ Orbitrap Velos™	25 000 to 100 000

Go to the next topic, "Selecting the Input Files for a New Analysis."

Selecting the Input Files for a New Analysis

Follow this procedure to select the input files for a new analysis. If you are reprocessing a previous analysis, the Files for Analysis area of the Analysis pane includes its input files. If you are starting a new analysis, the Files for Analysis area is empty.

* To select the input files that you want to process

- 1. Open the study of interest and start a new analysis.
- 2. On the Input Files page or the Samples page of the study, select the files of interest.
- 3. Right-click the selection and choose **Set As Input File**, or drag the files of interest to the Files for Analysis area of the Analysis pane (Figure 60).

Analysis Results Grouping & Ratios				Norkflo	Analysis 🗌 As Batch 🎲 Run 🛃 Save	e				
		Definiti			nput Files			Sample		
_	Samp 🔺		Sample Io	dentifier	Sample Type		Phenot	ype	Processing Step (Fully Processing) Edit	1
-				•	• •			•	Workflow:	
	S1	F1	blank		Blank	*	Lean	*	Result File: Enter result file name.	
	S2	F2	Pooled_do	dMS2	Identification	*	n/a	*	Resolutine. Enter resolutionen	
	S3	F3	ZDF_Fatty	1	Sample	*	Fatty	*	Files for Analysis: (0) 🛛 💥 Clea	ar Al
	S4	F4	ZDF_Fatty	_2	Sample	*	Fatty	•	La	
	S5	F5	ZDF_Fatty	_3	Sample	*	Fatty	*	Drag and the blank Input Files here	e
	S6	F6	ZDF_Lean	1	Sample	*	Lean	*	Colled_ddMS2	
	S7	F7	ZDF_Lean	_2	Sample	*	Lean	*	ZDF_Fatty_1	
	S8	F8	ZDF_Lean	_3	Sample	•	Lean	•	CDF_Fatty_2	
									CZDF_Fatty_3	
									CZDF_Lean_1	
									ZDF Lean 2	

Figure 60. Dragging input files to the Analysis pane

The file name of the last input file appears in the Result File box, and the Run command becomes available if the Workflows page contains a valid processing workflow. When you add more than one input file to the Files for Analysis area, the As Batch check box becomes available.

If the Caution symbol in the Processing Step title bar remains, the processing workflow contains an error.

Go to the next topic, "Naming the Result File for a New Analysis."

Naming the Result File for a New Analysis

After you add the files to be processed to the Analysis pane, the application automatically populates the Result File box with the file name of the first input file.

To change the name of the result file for an analysis

In the Analysis pane, select the default file name in the Result File box and type the new name.

Go to the next topic, "Setting Up the Sample Groups and Ratios for a New Analysis."

Setting Up the Sample Groups and Ratios for a New Analysis

Note When the processing workflow includes the Differential Analysis node, the Analysis Validation confirmation box opens if the analysis does not include defined sample groups and ratios on the Grouping & Ratios page.

***** To set up sample groups and ratios for a new analysis

- 1. Add input files to the Analysis pane (see "Selecting the Input Files for a New Analysis" on page 118).
- 2. Follow the instructions for the Sample Groups and Ratios page of the New Study and Analysis Wizard ("Setting Up the Sample Groups and Ratios for a New Analysis" on page 92).

Go to the next topic, "Troubleshooting Common Analysis Errors."

Troubleshooting Common Analysis Errors

After you set up an analysis, if a caution symbol appears in the Analysis pane and the Run button is unavailable, you must troubleshoot the analysis.

Tip The application does not know the polarity of the scan data or whether the data contains data-dependent or data-independent fragmentation (acquisition) scans. For best results, check the following:

- If the processing workflow contains any of these nodes—Create Mass Trace, Create FISh Trace, or Create Pattern Trace—verify that the node's ion polarity setting matches the data.
- If the processing workflow contains the Create FISh Trace node, verify that the setting for Fragmentation Mode matches the data.

✤ To troubleshoot an analysis

- 1. In the Analysis pane, point to the Caution symbol, 4, to display the list of missing analysis items, or check the error information in the Current Workflow Issues pane (Figure 62) below the Post-Processing pane on the Workflows page.
- 2. Using the information provided in the list of missing items, fix the analysis errors until the Caution symbol disappears (see "Common Analysis Errors" on page 122).

For example, Figure 61 shows the error message that appears before you customize the defined processing workflow provided in the Common Templates folder:

Degradants w Statistics Expected w FISh Scoring and Unknown ID w Online and Local Database Searches.cdProcessing WF

Analysis			🗆 As Batch 🞲 Run 📙 Save 🗙
Processing S	itep (Fully Processing)		Edit: 🔥
Workflow:	Degradants w Stats Expected w FISh Scoring	Create FISh Trace	Missing value for parameter 'Compound'
Result File:	Assay.cdResult	Search Mass Lists	Missing value for parameter 'Mass Lists'
▼ Files for	Analysis: (6)	Compound Class Scoring	Missing value for parameter 'Compound Classes
	-	Generate Expected Compounds	Missing value for parameter 'Compound'

Figure 61. Error messages for a defined processing workflow

Figure 62 shows the corresponding error messages in the Current Workflow Issues pane.

Figure 62. Current Workflow Issues pane

Current Workflow Issues					
Node Name	Issue Description	Parameter Name	Value		
Create FISh Trace	Missing value for parameter 'Compound'	Compound			
Search Mass Lists	Missing value for parameter 'Mass Lists'	Mass Lists			
Compound Class Scoring	Missing value for parameter 'Compound Classes'	Compound Classes			
Generate Expected Compounds	Missing value for parameter 'Compound'	Compound			

In addition, each workflow node that is missing a value for one of its parameters has an exclamation mark, ¹, in its upper right corner.

To fix the analysis errors for this example, you must select a compound for the Create FISh Trace node, one or more compounds for the Generate Expected Compounds node, and a mass list for the Search Mass Lists node.

Tip To view new compounds in the Compound list or new mass lists in the Input Files list, close the current analysis and start a new analysis. This means that you must select the processing workflow on the Workflows page and add the input files that you want to process from the Input Files page to the Analysis pane.

Go to "Submitting an Analysis to the Job Queue" on page 122.

Common Analysis Errors

To troubleshoot common analysis errors, see Table 19.

Note To troubleshoot validation issues, which occur after you submit a run, see "Common Validation Issues" on page 123.

Table 19. Common analysis errors

Error message	Error	Solution
The workflow does not contain a start node.	You have not added the Input Files node to the workflow.	Add the Input Files node to the beginning of the workflow.
No input files defined.	You have not added input files to the Files for Analysis area.	Add input files from the Input files or Samples page to the Files for Analysis area.
The current workflow does not contain any nodes.	You have not set up a processing workflow on the Workflows page.	Select or set up a processing workflow on the Workflows page.
<i>Node Name</i> Missing value for parameter " <i>Parameter Name</i> "	The processing workflow contains a node that requires a custom parameter setting.	Make the appropriate selections in the affected workflow nodes.
<i>Node Name</i> Missing connection for " <i>Connection Information</i> "	You have not connected node to the processing workflow.	Make the appropriate node connections.

Submitting an Analysis to the Job Queue

After you set up an analysis and troubleshoot any analysis errors, the Run command becomes available.

- To submit an analysis to the job queue
- 1. Decide whether you want to create one result file for the entire set of input files or one result file for each input file.
 - To create a single result file, leave the As Batch check box clear.
 - To create a separate result file for each input file, select the As Batch check box.
- 2. (Optional) In the Result File box, type a name for the result file to overwrite the default name.
- 3. In the Analysis command bar, click Run.

If the analysis contains no validation issues, the Job Queue page opens.

If the analysis contains issues, the Analysis Validation Issues prompt opens.

- 4. At the prompt, do the following:
 - a. Read the warning message.
 - b. Do one of the following:
 - To modify the analysis, click **Abort** and remedy the error.
 - To start data processing, click Ignore.

Related Topics

- Working with the Job Queue
- Common Validation Issues

Common Validation Issues

To troubleshoot common validation issues, see Table 20.

Table 20. Validation issues (Sheet 1 of 2)

Validation issue	Remedy	
The processing workflow includes the Differential Analysis node, but you have not set up sample groups and ratios.	• On the Sample Groups and Ratios page, set up the appropriate sample groups and ratios.	
	 On the Workflows page, delete the Differential Analysis node. 	
	-10-	
	• Do not change the analysis settings and click Ignore .	
The analysis includes sample groups and ratios, but the processing workflow does not include the Differential Analysis node.	On the Workflows page, add the Differential Analysis node to the workflow.	
IMPORTANT If the analysis does not includ node, the application does not run a differen Expected Compounds tables in the result file Ratios, Log2 Fold Change, P-value, and Adj	ntial analysis, and the Compounds and e does not include the following columns:	
The processing workflow includes Compound Identification or Pathway	On the Workflows page, add the Assign Compound Annotations node to the	
Mapping nodes, but does not include the Assign Compound Annotations node.	workflow.	
IMPORTANT If the analysis does not include the Assign Compound Annotations node, the application does not assign names or formulas to compounds in the Compounds table of a result file.		

 Table 20.
 Validation issues (Sheet 2 of 2)

Validation issue	Remedy
The processing workflow includes the Assign Compound Annotations node, but the processing workflow does not include any of the Compound Identification or Pathway Mapping nodes.	Click Ignore . –or– Add one or more Compound Identification
Mapping notes.	or Pathway Mapping nodes to the processing workflow.

Working with the Job Queue

When you submit an analysis for processing, the Job Queue page opens (Figure 63). You can also open the Job Queue page by choosing View > Show Job Queue from the menu bar.

The application can process two runs simultaneously. If you submit a second run while the first run is being processed, the status of the second run goes to Not Queued, Running, or Execution Failed. If you submit a third run while the application is processing the first two runs, its status goes to Waiting. If you pause a run that is waiting, its status goes to Sleeping.

Figure 63. Job queue with runs in various processing states

1	T Start Page × III Omeprazole Study × Z Job Queue ×					
ŵ	🏟 Pause 🏟 Resume 🎲 Abort 🐹 Remove 🥏 Refresh 👘 Open Results 🔲 Display Verbose Messages					
	Job Queue:					
	Execution State	Details	Progress	Туре	Name	Submitted at
- m						=
÷	Waiting		0 %	Processing	Assay 5	6/5/2018 4:47 PM
÷	Running		15 %	Processing	Assay 4	6/5/2018 4:40 PM
	Running		15 %	Processing	Assay 3	6/5/2018 4:40 PM
÷	Completed	Warnings	100 %	Processing	Assay 2	6/5/2018 4:39 PM

To work with the Job Queue page, see the following topics:

- Opening the Job Queue Page
- Pausing a Run That Is Waiting to be Processed
- Resuming a Paused Run
- Canceling a Run
- Removing Completed or Failed Runs from the Job Queue
- Refreshing the Job Queue
- Opening a Result File from the Job Queue
- Displaying Verbose Messages in the Job Queue

- Viewing the Processing Steps for a Job
- Filtering the Job Queue

Opening the Job Queue Page

The Job Queue page automatically opens when you submit a job for processing.

To open a hidden Job Queue page

From the Compound Discoverer window, choose **View > Show Job Queue**. The Job Queue page opens as a tabbed document.

Pausing a Run That Is Waiting to be Processed

You can only pause runs with a status of Waiting.

To pause a run that is waiting to be processed

- 1. In the Job Queue list, select the run that you want to pause.
- 2. In the command bar, click **Pause**.

The status of the run changes to Sleeping.

Resuming a Paused Run

✤ To resume a paused run

- 1. In the Job Queue list, select the paused run.
- 2. In the command bar, click **Resume**.

Canceling a Run

To cancel a run that is being processed

- 1. In the Job Queue list, select the run that you want to cancel.
- 2. In the command bar, click **Abort**.

When you cancel a run, its status changes to Aborted, and the application does not create a result file for the run.

Removing Completed or Failed Runs from the Job Queue

* To remove a completed or failed run from the Job Queue list

- 1. In the Job Queue list, select the run that you want to remove.
- 2. In the command bar, click **Remove**.

The selected run disappears from the Job Queue list. The application does not remove the result file from the study.

Refreshing the Job Queue

✤ To refresh the Job Queue list

In the command bar, click **Refresh**.

Opening a Result File from the Job Queue

* To open a result file from the Job Queue list

In the Job Queue list, do one of the following:

• Select the completed run of interest and click **Open Results** in the command bar.

-or-

• Double-click a completed run of interest.

The Results page for the selected run opens as a tabbed document in the application window.

Displaying Verbose Messages in the Job Queue

By default, the application only displays high-level processing information on the Job Queue page. To display all the available processing information, you can display the verbose messages.

* To display verbose messages in the Job Queue list

Select the **Display Verbose Messages** check box.

Viewing the Processing Steps for a Job

✤ To view the processing steps for a job

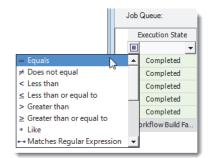
Click the expand icon, I, to the left of the job row.

Filtering the Job Queue

✤ To filter the Job Queue list

Note Use the filters for the column that you want to sort by. For example, to display only the runs that you ended before completion, follow this procedure.

1. Click the icon, 🗐, to the left of the Execution State filter list and select **Equals**.



2. In the Execution State filter list, select Aborted.

Job Queue:				
Execution State				
	=		•	
÷		(Custom)		
+		(Blanks)		
	_	(NonBlanks)		
- -	_	Aborted		
1	_	Completed		
+		Aborted		

The Job Queue list displays the canceled runs only.

3. To undo filtering, close and reopen the Job Queue page.

Job Queue Page Parameters

Table 21 describes the command bar and progress table on the Job Queue page.

 Table 21.
 Job Queue page features (Sheet 1 of 3)

Feature	Description
Command bar	
Pause	Pauses a job that is waiting to start.
	Selecting a job that is waiting to start activates this button.
Resume	Resumes a processing job that is sleeping.
	Selecting a job with a status of Sleeping activates this button.

Fastura	Description		
Feature	Description		
Abort	Stops processing and removes the selected job from the queue.		
	Selecting a job that is being processed activates this button.		
Remove	Selecting a completed job activates this button.		
Refresh	Refreshes the job queue list.		
Open Results	Opens the result file for the selected job.		
Display Verbose Messages	Displays more messages of less importance. When this check box is clear, the Job Queue displays no more than a few messages for each workflow node.		
Table columns			
Execution State	Displays the status of the job.		
	• Not Queued—The application takes a finite length of time to start a job after you click the Run command.		
	• Running—The application is currently processing the job. The application can process two jobs simultaneously. When you submit an analysis as a batch, each input file is processed as a separate job.		
	 Aborted—You canceled the job while the application was processing the analysis. 		
	• Execution Failed—The application was unable to complete the job.		
	• Waiting—The application has not begun to process the job.		
	 Sleeping—You paused the job while it was waiting in the queue. 		
	• Completed—The application has completed the analysis and you can open the result file.		
Details	Displays whether the job ran with or without warnings.		
Progress	Displays the progress of the run as a percentage.		
Туре	Displays the job type.		
Name	Displays the name of the result file.		
Submitted At	Displays the date and time when you submitted the run to the job queue.		
Study	Displays the name of the study.		

Table 21. Job Queue page features (Sheet 2 of 3)

Feature	Description
Data Source	Displays the location and file names of the input files for the current job.
Description	Displays the description that you typed in the Description box on the Workflows page.

Table 21. Job Queue page features (Sheet 3 of 3)

Reviewing or Reprocessing an Analysis

When you submit a run to the job queue, the run appears in the list of analyses on the Analysis Results page of the current study.

Use the Analysis Results page of a study to review or reprocess a completed analysis or to open a result file. An analysis consists of a processing workflow, optional sample groups and ratios, and the selected input files.

Note The Execution State on the Job Queue page updates more quickly than the Execution State on the Analysis Results page.

See these topics:

- Analysis Results Page Parameters
- Reviewing an Analysis
- Reprocessing an Analysis

Analysis Results Page Parameters

Table 22 describes the command bar and progress table on a study's Analysis Results page.

Table 22. Analysis Results page parameters (Sheet 1 of 2)

Feature	Description
Command bar	

Selecting an analysis in the list on this page activates these commands.

Open Result	Opens the selected result file.
Reprocess	Opens the Analysis pane with the list of input files that were used for the selected analysis. The Workflows page contains the processing workflow and the Grouping & Ratios page contains the sample groups and ratios used for the selected analysis.
Table columns	
Execution State	Displays the status of the analysis.

Feature	Description			
Creation Date	Displays the date and time when you submitted the run to the jo queue.			
File Name	Displays the file name of the result file.			
Description	Displays the description of the processing workflow that you or the originator typed in the Description box on the Workflows page.			
Below the table				
Show Associated Analysis/Hide Associated Analysis	Opens a locked Processing Step area that shows the names of the processing workflow and result file and a list of the input files for the analysis.			
Shortcut menu command	S			
Copy with Headers, Copy, Clear Selection, Cell Selection Mode, and Enable Row Grouping	See "Copying Table Entries to the Clipboard" on page 586 and "Grouping Table Rows" on page 582.			
Open Result	Opens the selected result file.			
Open Containing Folder	Opens the folder that contains the result file.			
Show Details	Opens the Analysis Sequence Details window where you can view the analysis processing workflow in the Workflow Tree pane and the Processing Step information in the Analysis pane. You cannot start runs from this window.			
Reprocess	Same functionality as the Reprocess command in the command bar.			

Table 22. Analysis Results page parameters (Sheet 2 of 2)

Reviewing an Analysis

Use the Analysis Sequence Details window to review an analysis.

✤ To review an analysis

1. On the Analysis Results page of a study, right-click the analysis and choose **Show Details**.

The Analysis Sequence Details window opens and displays the processing workflow in the Workflow Tree pane, and the result file name and input files in the Processing Step pane.

2. To view a workflow node's parameter settings, select it in the Workflow Tree pane to display its parameter settings in the Parameters pane on the left.

Note The Processing Step is locked, preventing you from reprocessing the analysis from this window.

Figure 64. Analysis Sequence Details window showing the parameter settings for the Detect Compounds node

🙈 Analysis Sequence Details		
Parameters of 'Detect Compounds'	👫 Open 🏢 Open Common 🛔 Save 📳 Save Common 🂝	Processing Step (Locked)
Show Advanced Parameters	Workflow: Untargeted Metabolomics with Statistics Detect Unkno	
1. General Settings Mass Tolerance [p. 5 ppm Intensity Tolerance 30 S/N Threshold 3 Min. Peak Intensity 1000000	Description: Untargeted Metabolomics workflow: Find and identify the differences between samples	Workflow: Untargeted Metabolomics with Statistics Detect Unknowns with ID using Online Databases Result File: C:\Tutorials\Metabolomics Study\ZDF Study\ZDF.cdResult
Ions [2M+ACN+H]+1; [2M+ACN+Na		▼ Files for Analysis: (8)
Min. Element Cour C H Max. Element Cou C90 H190 Br3 Cl4 K2 N10 Na2 O1	Align Retention Times	F1 blank Sample Type: [Blank], Phenotype: F2 Pooled_ddMS2 Sample Type: [Identification Only] F3 ZDF_Fatty_1 Sample Type: [Sample], Phenotype F4 ZDF_Fatty_2 Sample Type: [Sample], Phenotype F5 ZDF_Fatty_3 Sample Type: [Sample], Phenotype F6 ZDF_Fatty_3 Sample Type: [Sample], Phenotype F7 ZDF_Lean_1 Sample Type: [Sample], Phenotype F8 ZDF_Lean_3 Sample Type: [Sample], Phenotype III IV
Mass Tolerance [ppm] This parameter specifies the mass tolerance to be used for extracted ion chromatogram creation. Minimum value = 1 ppm Maximum value = 20 ppm	Detect Compounds Outrophysical Analysis Outrophysical Analysis The second	

Reprocessing an Analysis

You can reprocess an entire analysis or only the post-compound detection portions of an analysis.

To reprocess an analysis

- 1. On the Analysis Results page of a study, right-click the analysis and choose Reprocess.
 - If the Analysis pane is open and contains information for an analysis that you have not yet submitted to the job queue, the following prompt appears:

The analysis was not started. Do you really want to discard the analysis?

- If the Analysis pane is closed, it reopens with the settings from the selected analysis. In addition, the Grouping & Ratios tab and Workflows tab appear to the right of the study tabs.
- 2. If prompted, do the following:
 - a. Check the Analysis pane and decide whether you want to discard the current settings.

- b. Click **Yes** to replace the current settings in the Analysis pane with the settings from the selected analysis. Otherwise, click **No** to return to the in-progress analysis.
- 3. Do one of the following:
 - To reprocess the entire analysis, go to step 5.
 - To reprocess only part of the processing workflow, go to step 4.

Note You can reprocess any of the Peak Area Refinement nodes, Pathway Mapping nodes, Search nodes, and Compound Scoring nodes without reprocessing the entire workflow.

You must reprocess the entire workflow to reprocess any of the nodes for Spectrum Processing, Trace Creation, Compound Detection (except the Mark Background Compounds node), or Expected Compounds (except the Mark Background Compounds and Merge Features nodes).

- 4. To reprocess only part of the processing workflow, do the following:
 - a. Click the **Workflows** tab.

The processing workflow appears in the Workflow Tree (Reprocess) area. All the nodes are white with a gray tab, which indicates that they are not set for reprocessing.

b. Right-click the nodes that you want to reprocess and choose Reprocess.

The selected nodes and any related nodes revert to their original color, and the Run command becomes available in the Analysis pane.

Note If you are only partially reprocessing the existing input files, the Source File row appears in the analysis pane. You cannot edit the source file name.

c. If you are partially reprocessing the data, consider changing the result file name.

If you do not change the result file name, the application automatically appends a number to the file name when you submit the run.

- d. Go to step 7.
- 5. To reprocess the entire processing workflow, right-click the Input Files node and choose **Reprocess**. Then, at the prompt, click **OK**.
- 6. If necessary, make changes to the sample groups and ratios on the Grouping & Ratios page and the settings in the Analysis pane.

Note When you change any of the settings on the Grouping & Ratios page or the files for analysis in the Analysis pane, the application automatically reprocesses the entire processing workflow.

7. To start the analysis, click **Run**.

Analysis Pane Parameters

The Analysis pane appears to the right of the analysis pages when you start a new analysis or open an existing analysis template. Table 23 describes the parameters in the Analysis pane.

Table 23. Analysis pane parameters (Sheet 1 of 2)

Parameters	Description
Title bar	
As Batch check box	Available when you add more than one input file to the Files for Analysis area.
	• Clear (default setting)—The application creates one result file as it processes the input files in the Files for Analysis area.
	• Selected—The application creates one result file for each input file in the Files for Analysis area.
Run command	Submits the analysis to the job queue.
	Available when the Workflow Tree pane on the Workflows page contains a valid processing workflow and the Files for Analysis area contains a list of input files.
Save command	Opens the Save Analysis Template dialog box where you can provide a file name for the analysis template and save it to an appropriate directory.
×	Closes the Analysis pane, the Workflows page, and the Grouping & Ratios page.
Processing Step area	
Edit	Opens the Workflows page.
<u>^</u>	If the analysis includes errors, such as missing parameter settings or no input files, a Caution symbol appears to the far right of the Processing Step. To display the error list, point to the Caution symbol.
Workflow	By default, the text matches the text in the Workflow box on the Workflows page. You cannot change the text in the Analysis pane To change the name of the processing workflow, edit the text on the Workflows page.
Result File	Specifies the file name for the result file. This box is empty until the Files for Analysis area lists at least one input file.
	The name of the first input file automatically populates the Resul File box. If the analysis creates only one result file, you can type a name for the result file in the Result File box.

Parameters	Description
Source File	Displays the filename of the original result file.
	Available when you set up an analysis for partial reprocessing.
Files for Analysis: (#)	Displays the number of input files in the Files for Analysis area.
Files for Analysis area	Displays the names of the input files after you add the input files from the Input Files page or the Samples page of the study to this area.

Table 23. Analysis pane parameters (Sheet 2 of 2)

Related Topics

- Common Analysis Errors
- Selecting the Input Files for a New Analysis

Creating and Editing Processing Workflows

These topics describe how to create and edit processing workflows on the Workflows page of a study.

Contents

- Processing Workflows
- Workflows Page
- Editing a Processing Workflow
- Setting Up Isotope Patterns for Pattern Scoring or Pattern Traces
- Creating an Isotope Patterns List
- Extracting Analog Traces
- Extracting PDA Traces
- Creating a Completely New Processing Workflow
- Connecting the Workflow Nodes
- Input and Output Nodes for the Workflow Nodes
- Saving a Processing Workflow as a Template

 $\mathbf{\hat{h}}$

Processing Workflows

The Compound Discoverer application uses a processing workflow to analyze the MS data acquired with a high-resolution, accurate-mass LC/MS/MS instrument. In addition to analyzing the MS data, the application can display chromatograms acquired with a photo-diode array (PDA) or ultraviolet-visible (UV-Vis) detector that is controlled by a Thermo Scientific data system or an analog detector that is connected to the mass spectrometer's analog input channels.

Workflows Page

Within an existing study, use the Workflows page to create or edit processing workflows. The Workflows page is a tabbed page that opens to the right of the tabbed study pages when you start a new analysis or reprocess an existing analysis. You cannot edit a processing workflow from outside a study.

For information about opening the Workflows page and its command bar and shortcut menu, see these topics:

- Opening the Workflows Page
- Workflows Page Command Bar
- Workflows Page Shortcut Menu

Opening the Workflows Page

The Workflows page appears to the right of the four study pages and next to the Grouping and Ratios page after you set up a new study and a new analysis with the New Study and Analysis wizard.

Closing a study erases the analysis settings, and reopening a study only opens the study pages.

* To open the Workflows page

Do one of the following:

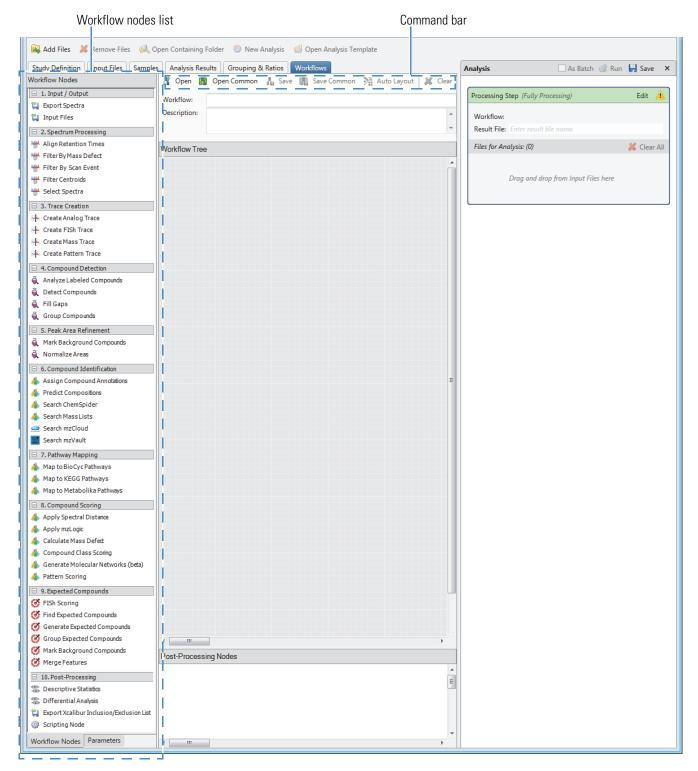
- Use the New Study and Analysis wizard to set up a new study. Then, click the **Workflows** tab.
- Open an existing study, click **New Analysis** in the Study command bar, and then click the **Workflows** tab.

-or-

• From the Analysis Results page of a study, select an analysis, click **Reprocess**, and then click the **Workflows** tab.

The Workflow Tree and Workflow Nodes panes appear (Figure 65). For a new analysis, the Workflow Tree pane is empty. For an existing analysis, the Workflow Tree pane usually contains a processing workflow.

Figure 65. Workflows page without a selected processing workflow



Workflows Page Command Bar

Table 24 describes the Workflows page commands.

Table 24. Workflows page commands

Command	Description
Open	Opens the Open Workflow dialog box for locating and opening a processing workflow.
Open Common	Opens the Open Workflow dialog box to the following folder where the application installs the three common processing workflow templates:
	C:\Users\Public\Public Documents\Thermo\Compound Discoverer 3.1\Common Templates
Save	Opens the Save Workflow dialog box for selecting a folder and entering a file name for the processing workflow in the Workflow Tree pane.
Save Common	Opens the Save Workflow dialog box to the folder where the application installs the common processing workflow templates. Saves the current processing workflow in the Workflow Tree pane to the Common Templates folder.
Auto Layout	Automatically formats the layout of the workflow nodes.
Clear	Clears the Workflow Tree pane.

Workflows Page Shortcut Menu

Table 25 describes the commands in the shortcut menu for the Workflow Tree pane or a workflow node.

Table 25. Workflow Tree pane and workflow node shortcut menu commands (Sheet 1 of 2)

Command	Description			
Cut	Removes the node from the workflow.			
	Selecting any workflow node enables this command.			
Copy/Paste	Adds a copy of the selected node to the Workflow Tree pane.			
	Selecting a workflow node that can appear more than once in a workflow enables this command.			
Auto Layout	Automatically formats the layout of the workflow nodes.			

Command	Description	
Hide Node Numbers	Turns the numbering on or off.	
Reprocess	Sets the selected node for reprocessing or sets the selected node and its related nodes for reprocessing.	
	Available for an analysis that you have set up for reprocessing [Workflow Tree (Reprocess)].	

Table 25. Workflow Tree pane and workflow node shortcut menu commands (Sheet 2 of 2)

Editing a Processing Workflow

You can modify a processing workflow by adding and deleting workflow nodes and by changing the parameter settings in the workflow nodes.

The application automatically connects some of the workflow nodes as you add them to a processing workflow. But for some of the workflow nodes, you must make the appropriate connections. When a node is missing a connection, a Caution symbol appears in its upper right corner.

Several workflow nodes require custom parameter selections. When a workflow node is missing a custom parameter selection, an exclamation mark appears in its upper right corner.

Follow these topics as needed.

- Fixing a Workflow Node That Has a Caution Symbol
- Fixing a Workflow Node that Has an Exclamation Mark
- Adding a Workflow Node to a Processing Workflow
- Deleting a Workflow Node from a Processing Workflow
- Editing the Parameter Settings for a Workflow Node

Fixing a Workflow Node That Has a Caution Symbol

When the node is missing a connection, a Caution symbol, \bigtriangleup , appears in its upper right corner. If the node is also missing a required parameter selection or the selected item is not found in its corresponding list or library, an exclamation mark appears after you fix the missing connection.

* To fix a workflow node that is labeled with a Caution symbol

1. To view the validation errors, point to the Caution symbol in the Analysis pane.

A missing connection begins with the following text:

Missing connection for

2. Make the appropriate node connections.

Related Topics

- Input and Output Nodes for the Workflow Nodes
- Connecting the Workflow Nodes

Fixing a Workflow Node that Has an Exclamation Mark

To edit a processing workflow, you must open it from the Workflows page within a study.

An exclamation mark, (), appears in the upper right corner of a node when the node is missing a required parameter selection or the selected item is not found in its corresponding list or library.

To fix a workflow node that is labeled with an exclamation mark

1. To view the validation errors, point to the Caution symbol in the Analysis pane or review the issues described in the Current Workflow Issues pane below the Post-Processing Nodes pane.

For a missing parameter selection, the issue description includes the following text:

Missing value for parameter 'Node'

- 2. To fix the workflow node error, select the node in the Workflow Tree pane.
- 3. In the Parameters of '*Workflow Node*' page to the left, make a selection for the missing parameter value.

Related Topics

Common Analysis Errors

Adding a Workflow Node to a Processing Workflow

To edit a processing workflow, you must open it from the Workflows page within a study.

To add a node to the processing workflow

1. Select the node in the Workflow Nodes pane and drag it to the Workflow Tree pane.

2. If necessary, make the appropriate node connections.

Note A Caution symbol appears in the upper right corner of a workflow node that is missing a connection. Because a missing connection takes priority over a missing parameter setting, if a workflow node is also missing a parameter setting, a blue circle with an exclamation mark appears after you fix the missing connection.

3. To display a node's parameters, click the node in the Workflow Tree pane.

The Parameters page lists the parameters for the selected node.

4. Edit the parameter settings for the node.

Related Topics

- Connecting the Workflow Nodes
- Input and Output Nodes for the Workflow Nodes

Deleting a Workflow Node from a Processing Workflow

To edit a processing workflow, you must open it from the Workflows page within a study.

- To delete a node from a processing workflow
- 1. Right-click the node in the Workflow Tree pane and choose Cut.
- 2. Check whether any validation issues appear.
- 3. (Optional) To save the processing workflow for reuse, click **Save**. Then, rename the workflow if necessary, select an appropriate folder, and click **Save**.

Tip If you use the processing workflow for the current analysis, you can reuse the processing workflow without saving it. After you run the analysis successfully, you can rerun the analysis by selecting the completed analysis on the Analysis Results page of a study.

Editing the Parameter Settings for a Workflow Node

To edit a processing workflow, you must open it from the Workflows page within a study.

To edit the parameter settings for a workflow node

1. In the Workflow Tree pane, select the node.

The Parameters page opens with the parameters for the selected node.

2. Click Show Advanced Parameters below the Parameters page title bar.

If the node contains hidden advanced parameters, the advanced parameters appear below the basic parameters.

When you place the cursor in the box to the right of the parameter name, information about the parameter appears at the bottom of the Parameters page.

3. Enter the appropriate values or make the appropriate selection for each parameter.

Related Topics

See Chapter 7, "Workflow Nodes."

Setting Up Isotope Patterns for Pattern Scoring or Pattern Traces

Use the Isotope Ratio Editor dialog box to set up the pattern and the required isotopes for the Create Pattern Trace node or the Pattern Scoring node.

You can access the Isotope Ratio Editor dialog box from the Create Pattern Trace node or the Pattern Scoring node.

To set up an isotope pattern, follow these topics as needed:

- Opening the Isotope Ratio Editor
- Defining an Isotope Pattern from an Elemental Composition
- Copying an Elemental Composition from the Expected Compounds Library
- Exporting a Mass Spectrum from a RAW File to the Clipboard
- Isotope Ratio Editor Parameters

Opening the Isotope Ratio Editor

To open the Isotope Ratio Editor

- 1. Open or create a processing workflow on the Workflows page.
- 2. Add one or both of these nodes to the processing workflow: Create Pattern Trace or Pattern Scoring.
- 3. Do one of the following:
 - a. Click the Create Pattern Trace node.

The parameters for the node appear in the Parameters pane.

- b. To display the browse icon, click the **Isotope Ratios** box.
- c. Click the browse icon,

The Isotope Ratio Editor dialog box opens (Figure 66).

-or-

a. Click the **Pattern Scoring** node.

The parameters for the node appear in the Parameters pane.

- b. To display the browse icon, click the **Isotope Ratios** box.
- c. Click the browse icon,

The Pattern List Editor dialog box opens.

d. In the Pattern Editor dialog box, click Add Patterns.

The Isotope Ratio Editor dialog box opens (Figure 66).

Figure 66. Isotope Ratio Editor dialog box with an isotope pattern for a defined elemental composition

Isotope Ratio	o Editor	×		
Isotope Ratio Definition Type				
O Define f	from Elemental Composition Formula			
C Custom	n Isotope Ratio Pattern			
Elemental Co	iomposition: C17 H19 N3 O3 S			
Int. Thre	reshold [%]: 0.10 Resolution: 60000 Charge: 1 📩			
150 - ⊮ ⊉ 100 -	345.11416			
0 - 000 - 0	346.11748 347.11001 348.11337 346.11748 347.12019 348.12248			
	345 346 347 348 m/z	349		
Denvind				
Required	Mass Shift Intensity [%]	100.00		
▼ ▼	1.0032	19.14		
	1.99585	4.57		
	2 00603			
		2.17		
	2.99920	0.87		
	3.00832	0.20		
	ОКС	ancel		

Defining an Isotope Pattern from an Elemental Composition

- * To set up the pattern for a compound by using its elemental composition
- 1. In the Isotope Ratio Definition Type area of the Isotope Ratio Editor dialog box, select the **Define from Elemental Composition Formula** option.
- 2. In the Elemental Composition box, type or paste the alphanumeric elemental composition of the compound of interest.

The application automatically populates the Mass Shift and Intensity [%] columns.

Tip To enter an elemental composition for a labeled compound, use brackets to identify the type and number of labeled atoms for each element.

For example, to enter the elemental composition of omeprazole where only one of the carbon atoms has been replaced with carbon-13, type C16 [13]C H19 N3 O3 S.

- 3. Do the following:
 - In the Int. Threshold [%] box, type the relative intensity threshold.

The application removes isotopic peaks below this relative intensity threshold from the Mass Shift versus Intensity [%] table.

• In the Resolution box, type the resolution for the scans.

Note The scan header in the raw data file lists the resolution of each scan, and the instrument method that is associated with the raw data file lists the resolution of each scan event.

• In the Charge box, type or select the charge state of the ions.

The application uses the charge state to display the theoretical mass spectrum in the graphical display.

4. Select the check boxes of the required peaks.

Copying an Elemental Composition from the Expected Compounds Library

- To copy a compound's elemental composition from the Expected Compounds library
- 1. If the Isotope Ratio Editor dialog box is open, close it.
- 2. From the menu bar, choose Libraries > Expected Compounds

The Expected Compounds library opens.

3. Right-click anywhere on the page and choose Cell Selection Mode.

4. Click the elemental composition of interest.

The table cell turns a lighter blue than the remaining cells in the table row.

- 5. Right-click the highlighted cell and choose **Copy**.
- 6. Open the Isotope Ratio Editor dialog box.
- 7. Right-click in the Elemental Composition box and choose Paste.

Setting Up a Custom Isotope Pattern

✤ To set up a custom pattern

1. In the Isotope Ratio Editor dialog box, select the **Custom Isotope Ratio Pattern** option.

Below the Isotope Ratio Definition Type area, the available parameters change (Figure 67).

Figure 67. Custom Isotope Ratio Pattern view

Isotope F	Ratio Edito	r					×
Isotop	Isotope Ratio Definition Type						
C De	fine from E	lemental (Composition	Formul	а		
⊙ Cu	stom Isotop	be Ratio F	attern				
Ge	t from com	position f	ormula		Get from (clipboard	
Intensity [%]		0.2	0.4	0.6 m/z	0.8	1.0	1.2
	Mass Shit	t	Intensi	ty [%]			
Þ	0		100				
*							
					ОК	Ca	ancel

- 2. To set up the custom isotope ratio pattern, do one of the following:
 - Type values in the Mass Shift and Intensity boxes.
 - Click Get from Composition Formula.

The application uses the text in the hidden Elemental Composition box.

• Click Get from Clipboard.

The application uses the mass spectrum data that you copied to the Clipboard.

Exporting a Mass Spectrum from a RAW File to the Clipboard

* To export a mass spectrum from a raw data file to the Clipboard

- 1. From the FreeStyle[™] data-visualization application, do the following:
 - a. Open the raw data file that contains the mass spectrum of interest and make the mass spectrum view or the spectrum list view the active view.
 - b. In the Exports area of the Spectrum Workspace Options toolbar, click Selection AS.
 - c. In the Copy to Clipboard/Export dialog box, select the **To CSV File** option and click **OK**.
 - d. In the Export Data dialog box, select a folder, name the file, and click Save.

The spreadsheet opens in a spreadsheet application.

2. In the spreadsheet application, select up to 20 rows of m/z and intensity values and copy them to the Clipboard. Do not select the spectrum header information.

Isotope Ratio Editor Parameters

Table 26 describes the parameters in the Isotope Ratio Editor dialog box.

Table 26. Isotope Ratio Editor dialog box parameters (Sheet 1 of 3)

Parameter	Description		
Isotope Ratio Definition Type			
These two options defin Type area.	e the parameters that appear below the Isotope Ratio Definition		
Define from Elemental	Selecting this option makes the following features visible:		
Composition Formula	Elemental Composition box		
	• Int. Threshold [%] box		
	• Resolution box		
	Charge box		

Parameter	Description			
Custom Isotope Ratio	Selecting this option makes the following features visible:			
Pattern	Get from Composition Formula button			
	Get from Clipboard button			
	• Bar graph of Intensity [%] versus the <i>m/z</i> ratio of the isotopic mass peaks			
	• Data entry table where you define a custom isotope pattern in terms of the mass shift and intensity of each mass peak			
	Use this option to create a custom isotope pattern.			

 Table 26.
 Isotope Ratio Editor dialog box parameters (Sheet 2 of 3)

Elemental composition view

Selecting the Define from Elemental Composition Formula option makes the following parameters visible: Elemental Composition, Int. Threshold [%], Resolution, and Charge.

Elemental Composition	Specifies the elemental composition of the compound of interest.
	When you type a composition in this box, the application automatically creates a table of mass shifts and intensities.
	Default: Empty (unless you have already specified the elemental composition in the Isotope Ratios box under General Settings for the Pattern node)
Int. Threshold [%]	Specifies the intensity threshold of the isotope pattern.
	Default: 0.10; range: 0–100
Resolution	Specifies the resolution of the isotope pattern.
	Default: 60 000; range: 2–1 000 000 000
Charge	Specifies the charge of the ion fragment.
	Default: 1; range: 1–100

Graph of Intensity [%] versus m/z value

Displays a graph of the full isotope distribution in the mass shift and intensity table.

Table 26. Isotope Ratio Editor dialog box parameters (Sheet 3 of 3)

Parameter	Description	
Mass shift and inte	nsity table	

Use this table to specify the required peaks in the isotope pattern.

When you select the Define from Elemental Composition Formula option, the application automatically populates this table; you cannot edit the entries.

Required (for an isotope pattern defined from a user-specified elemental composition)	Specifies whether the isotope is required.When the check box is selected, the isotope is required.When the check box is clear, the isotope is not required.
Mass Shift	Specifies the mass shift from the pattern base peak (A0).
Intensity [%]	Specifies the relative intensity [%] of the isotope to the pattern base peak.
Rows	Specify the values for the isotopes.

Custom isotope ratio pattern view

Selecting the Custom Isotope Ratio Pattern option makes the following buttons visible: Get from Composition Formula and Get from Clipboard.

Get from Composition Formula	Creates an isotope pattern for an elemental composition formula. The application reads the elemental composition that you entered in the Elemental Composition box before selecting the Custom Isotope Ratio Pattern option.
Get from Clipboard	Imports the isotope pattern from the Clipboard. You can export a custom pattern to the Clipboard from the spectrum list view in the Qual Browser application or from a third-party software application.

Graph of Intensity [%] versus m/z value

Displays a graph of the full isotope distribution in the mass shift and intensity table.

Mass shift and intensity table

When you select the Custom Isotope Ratio Pattern option, you can edit the mass shift and intensity values for the isotope pattern.

Creating an Isotope Patterns List

The Pattern Scoring node requires a list of isotope patterns.

* To add isotope patterns to the Isotope Patterns node

Do either of the following:

- a. In the node's Isotope Patterns box, type the elemental composition formulas for the defined isotope patterns that you want to compare. Separate each pattern with a semicolon and a space—for example, C15S; C17S.
- b. Click the browse icon,, to the right of the Isotope Patterns box of the Isotope Patterns node.

The Pattern List Editor opens (Figure 68).

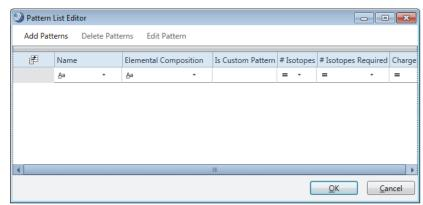
- c. Select the isotope pattern that you want to edit, and then click Edit Pattern.
- d. To edit the pattern, see "Setting Up Isotope Patterns for Pattern Scoring or Pattern Traces" on page 142.

-or-

a. Click the browse icon,, to the right of the Isotope Patterns box of the Isotope Patterns node.

The Pattern List Editor opens (Figure 68).

Figure 68. Pattern List Editor



b. For each pattern that you want to add, click **Add Patterns**, and then set up the isotope pattern in the Isotope Ratio Editor.

After you add a pattern, it appears in the Patterns table of the Pattern List Editor dialog box. The Name and Elemental Composition columns display the elemental composition of the pattern. The text in the Name column is editable.

c. Click **OK** to close the Pattern List Editor dialog box.

The isotope patterns appear in the Isotope Patterns box of the Isotope Patterns node.

Table 27 describes the commands and table columns in the Pattern List Editor dialog box.

Parameter	Description
Commands	
Add Patterns	Opens the Isotope Ratio Editor dialog box for setting up an isotope pattern.
Delete Patterns	Deletes the selected pattern.
Edit Pattern	Opens the Isotope Ratio Editor dialog box for editing the selected pattern.
Columns	
Name	By default, displays either the elemental composition for each isotope pattern that you create with a defined elemental composition or the text " <i>CustomPattern #</i> " for each isotope pattern that you create by setting up a custom isotope ratio pattern.
	You can edit the text in this column.
Elemental Composition	Displays the elemental composition for each isotope pattern that is based on a defined elemental composition.
Is Custom Pattern	Displays an X for each isotope pattern that is based on a custom isotope ratio pattern.
# Isotopes	Displays the number of isotopes in the defined pattern that are above the intensity threshold.
# Isotopes Required	Displays the number of isotopes required to calculate the SFit% score.

 Table 27.
 Pattern List Editor parameters

Related Topics

• Setting Up Isotope Patterns for Pattern Scoring or Pattern Traces

Extracting Analog Traces

To extract analog traces from your raw data, add the Create Analog Trace node to the processing workflow. Analog traces include data from a UV-Vis detector, analog data from LC devices controlled by the Xcalibur data system or an equivalent Thermo Scientific application, and analog signals from devices connected to the analog channels on the communications panel of your mass spectrometer.

To extract UV or analog traces from the raw data

- 1. Under General Settings, select UV or Analog from the Trace Type list.
- 2. In the RT offset [min] box, type the offset time, in minutes, for the UV-Vis or Analog trace.

If there is a time difference between when the sample enters the mass spectrometer and the UV-Vis or analog detector's flow cell, use the offset time to align the chromatographic traces. A value of 0 indicates that the UV-Vis or analog detector and the mass spectrometer detected the sample simultaneously.

3. In the Custom Label box, type text to identify the trace in the Specialized Traces table of the result file window.

Related Topics

• Create Analog Trace Node

Extracting PDA Traces

To extract the traces acquired by a photodiode array detector from your raw data, add the Create Analog Trace node to the processing workflow.

- To extract a PDA trace from the raw data
- 1. In the General Settings area, do the following:
 - In the Trace Type list, select **PDA**.
 - In the RT offset [min] box, type the offset time, in minutes, for the PDA traces.

If there is a time difference between when the sample enters the mass spectrometer and the PDA detector's flow cell, use the offset time to align the chromatographic traces. A value of 0 indicates that the PDA detector and mass spectrometer simultaneously detect the sample.

• In the Custom Label box, type text to identify the trace in the Specialized Traces table on the result file page.

- 2. In the PDA Settings area, do the following:
 - To extract a plot of the average intensity of the scanned wavelength range versus time, select **True** for Total Scan.

This trace is labeled as a Total Scan in the Specialized Traces table.

• To extract a plot of the spectrum maximum of the scanned wavelength range versus time, select **True** for Spectrum Maximum.

This trace is labeled as a Spectrum Maximum in the Specialized Traces table.

• To extract a plot for a specified wavelength range, select **True** for Wavelength Range. Then, type the wavelength range in the Min. and Max. Wavelength boxes.

Related Topics

• Create Analog Trace Node

Creating a Completely New Processing Workflow

A processing workflow is part of an analysis, and you can only perform analyses from inside a study. Therefore, to edit or create a processing workflow, you must open a study and start a new analysis or open an analysis template.

A processing workflow always begins with the Input Files node. All processing workflows that process MS data require the Select Spectra node. The only processing workflow that does not require the Select Spectra node is limited to processing the data from an analog detector.

- To create a completely new processing workflow from an empty Workflows page
- 1. Open the Workflows page. If the Workflows Tree area contains a processing workflow that you do not want to edit, click **Clear**.
- 2. Drag the required **Input Files** node from the Workflow nodes pane to the Workflow Tree pane.

The Input Files node reads the information in the raw data files.

3. To process the spectral data, drag the Select Spectra node to the Workflow Tree pane.

The Input Files node automatically connects to the Select Spectra node. The Select Spectra node reads and filters the MS scan data in the raw data files. The default parameter settings pass all the scan data to the next node.

4. To align multiple input files, drag the **Align Retention Times** node to the Workflow Tree pane. Then, connect the Select Spectra node to it.

The Align Retention Times node chromatographically aligns the MS scan data in the input files.

- 5. To find expected compounds, do the following:
 - a. Check whether your Expected Compounds library contains the compounds of interest (see "Modifying the Expected Compounds List" on page 498).
 - b. Drag the **Find Expected Compounds** node to the Workflow Tree pane. Then, connect the Align Retention Times node to it.

Note The Find Expected Compounds node accepts input from any of the Spectrum Processing nodes.

IMPORTANT For analyses with multiple input files, always add the Align Retention Times node to the processing workflow.

- c. Drag one or more **Generate Expected Compounds** nodes to the Workflow Tree pane.
 - To apply different transformation rules to multiple compounds, drag multiple Generate Expected Compounds nodes to the Workflow Tree pane, one for each set of rules.
 - To apply the same transformation rules to one or more compounds, drag a single Generate Expected Compounds node to the Workflow Tree pane.

IMPORTANT The Generate Expected Compounds node generates a list of expected compounds by using one or more user-specified library compounds and a set of user-specified chemical reactions.

The Compounds parameter is empty until you select the compounds of interest. If you submit an analysis to the job queue without selecting the compounds of interest, a Caution symbol appears.

- d. For each Generate Expected Compounds node, select the compounds of interest (see "Generate Expected Compounds Node" on page 234).
- e. Connect the Generate Expected Compounds node or nodes to the Find Expected Compounds node.
- f. Drag the Group Expected Compounds node to the Workflow Tree pane.

The Find Expected Compounds node automatically connects to it.

g. To add FISh scoring to the processing workflow for targeted compounds, drag the **FISh Scoring** node to the Workflow Tree pane.

The Group Expected Compounds node automatically connects to it.

Note FISh scoring adds a significant amount of processing time to an analysis.

- h. Do one of the following:
 - To detect unknown compounds in addition to the expected compounds, go to step 6.
 - To merge the features found by the Expected Compounds node and detected by the Unknown Compounds node, go to step 10.
- 6. To detect unknown compounds, do the following:
 - a. Drag the **Detect Compounds** node to the Workflow Tree pane. Then, connect the Align Retention Times node to the Detect Compounds node.
 - b. Drag the Group Compounds node to the Workflow Tree pane.

The Detect Compounds node automatically connects to it.

7. To identify unknown compounds, do the following:

Note The Group Compounds node automatically connects to all the Compound Identification nodes. In addition, the Group Expected Compounds node automatically connects to the Search Mass Lists, Search mzCloud, and Search mzVault nodes.

a. Drag the **Predict Compositions** node to the Workflow Tree pane.

IMPORTANT For best results, always include the Predict Compositions node in a processing workflow for untargeted compounds. Without the Predict Compositions node, the workflow does not report the elemental compositions of the unknown compounds without a hit from the online search databases or the local Metabolika database.

- b. To assign a name, formula, and structure to the compounds in the Compounds table, drag the **Assign Compound Annotations** node to the Workflow Tree pane.
- c. (Optional) To search the online ChemSpider database, drag the **Search ChemSpider** node to the Workflow Tree pane.
- d. (Optional) To search mass lists, drag the **Search Mass Lists** node to the Workflow Tree pane. In the Workflow Tree pane, select the **Search Mass Lists** node, and then, in the parameters pane, select the appropriate mass lists (see).
- e. (Optional) To search the online mzCloud database, drag the **Search mzCloud** node to the Workflow Tree pane.
- f. (Optional) To search your local mzVault database, drag the **Search mzVault** node to the Workflow Tree pane. In the Workflow Tree pane, select the **Search mzVault** node, and then, in the parameters pane, select libraries from the mzVault Library list.

8. To map detected compounds to a biochemical pathway, do any of the following:

Note The Group Compounds node automatically connects to all the Pathway Mapping nodes.

- To search the local Metabolika pathways, drag the **Map to Metabolika Pathways** node to the Workflow Tree pane.
- To search the KEGG database, drag the **Map to KEGG Pathways** node to the Workflow Tree pane. The Group Compounds node automatically connects to it.
- To search the BioCyc database, drag the **Map to BioCyc Pathways** node to the Workflow Tree pane.

The Group Compounds node automatically connects to it.

Note For information about adding the Mark Background Compounds nodes to an untargeted workflow, see "Marking Background Compounds in an Untargeted Analysis" on page 28. For information about using the Normalize Areas node for batch normalization, see "Quality Control Samples for Batch Normalization" on page 34.

 To rank the hits from the Search ChemSpider node, Map to Metabolika Pathways node, Search Mass Lists node, and Map to BioCyc Pathways node, drag the Apply mzLogic node to the Workflow Tree pane.

When present, the Search ChemSpider node, Map to Metabolika Pathways node, Search Mass Lists node, and Map to BioCyc Pathways node automatically connect to it.

 To compare the features found by the Find Expected Compounds and Detect Compounds nodes if applicable, drag the Merge Features node to the Workflow Tree pane.

The Find Expected Compounds and Detect Compounds nodes automatically connect to it.

- 11. To add Trace Creation nodes to the processing workflow, do the following:
 - For each UV, PDA, or analog trace that you want to extract, drag a **Create Analog Trace** node from the Workflow Nodes pane to the Workflow Tree pane.

The Input Files node automatically connects to it.

Note In this example, you connect the Align Retention Times node to the Create Pattern Trace, Create FISh Trace, and Create Mass Trace nodes. These nodes accept input from any of the Spectrum Processing nodes.

• For each pattern trace that you want to extract, drag a **Create Pattern Trace** node to the Workflow Tree pane. Then, connect the Align Retention Times node to each Create Pattern Trace node.

- For each Fragment Ion Search (FISh) trace that you want to create, drag a **Create FISh Trace** node to the Workflow Tree pane. Then, connect the Align Retention Times node to each Create FISh Trace node.
- For each mass trace that you want to extract, drag a **Create Mass Trace** node to the Workflow Tree pane. Then, connect the Align Retention Times node to each Create Mass Trace node.
- 12. In the Post-Processing Nodes pane, do the following:
 - To add a differential analysis (volcano plot) to the workflow for the specified sample groups and ratios, drag the **Differential Analysis** node to the Post-Processing Nodes pane below the Workflow Tree pane.
 - To add the descriptive statistics columns to the Compounds and Expected Compounds tables in the result file, drag the **Descriptive Statistics** node to the Post-Processing Nodes pane below the Workflow Tree pane. The descriptive statistics columns provide information about the chromatographic peak areas (mean area, median area, minimum area, and so on) for each detected compound.

Note In a result file, the descriptive statistics columns are hidden by default.

- 13. To export the MS scan data in the raw data files to a common data format, drag an **Export Spectra** node to the Workflow Tree pane. Then, connect one of these nodes to the Export Spectra node:
 - Select Spectra
 - Align Retention Times

Note The Export Spectra node does not export analog data.

Related Topics

- Workflow Nodes
- Connecting the Workflow Nodes

Connecting the Workflow Nodes

Although most of the processing workflow nodes automatically connect to the appropriate input and output nodes, you might need to manually connect some of the nodes.

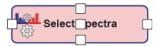
For a custom processing workflow, you must manually connect these nodes if you add them to the workflow:

- Spectrum Processing nodes
- Find Expected Compounds node
- Trace Creation nodes (except for the Create Analog Trace node)

To connect the workflow nodes

1. Point to the input node of interest.

Five white boxes appear, with one box at the center of the node and the other boxes at the center of each side.



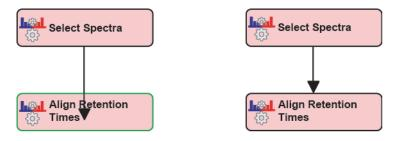
2. Click one of the white boxes and hold down the mouse button until a red border and an arrowhead appear.



3. Continue holding down the mouse button as you drag the arrowhead to the output node of interest.

Depending on the compatibility, one of the following occurs:

- If the selected input node is not compatible with the output node, a red border appears around the output node.
- If the selected input node is compatible with the output node, a green border appears around the output node. When you release the mouse button, a directional arrow connects the input node to the output node.



Note The Input Files node automatically connects to the Create Analog Trace node and the Select Spectra node.

4. To automatically format the layout of the workflow nodes, click **Auto Layout** in the Workflows command bar.

You can use the following nodes multiple times in a processing workflow: the Generate Expected Compounds node, all the Trace Creation nodes, and all the Spectrum Processing nodes, except for the Align Retention Times node.

Related Topics

- Workflows Page Command Bar
- Input and Output Nodes for the Workflow Nodes

Input and Output Nodes for the Workflow Nodes

Table 28 describes the node connections.

 Table 28. Input and output nodes for each workflow node (Sheet 1 of 3)

Workflow node	Input nodes	Output nodes
1.0 Input/Output		
Export Spectra Node	Select Spectra	Filter Centroids
Input Files (Begins every processing workflow.)	None	Select SpectraCreate Analog Trace
2.0 Spectrum Processing		
Align Retention Times Node	Spectrum Processing	 Spectrum Processing Trace Creation Find Expected Compounds Detect Compounds
Filter Centroids Node	Spectrum Processing	 Spectrum Processing Trace Creation Find Expected Compounds Detect Compounds
Filter By Mass Defect Node	Spectrum ProcessingGenerate Expected Compounds	 Spectrum Processing Trace Creation Find Expected Compounds Detect Compounds
Filter By Scan Event Node	Spectrum Processing	 Spectrum Processing Trace Creation Find Expected Compounds Detect Compounds
Select Spectra Node	Input Files	 Spectrum Processing Trace Creation Find Expected Compounds Detect Compounds
3.0 Trace creation		
Create Analog Trace Node	Input Files	None
Create Mass Trace Node	Spectrum Processing	None
Create FISh Trace Node	Spectrum Processing	None
Create Pattern Trace Node	Spectrum Processing	None
4.0 Compound Detection		
Analyze Labeled Compounds Node	Assign Compound Annotations	None

Detect Compounds NodeSpectrum Processing· Group Compounds · Merge FeaturesFill Gaps NodeGroup Compounds· Normalize Areas · Mark Background CompoundsGroup Compounds NodeDetect Compounds· Compound Identification · Pathway Mapping · Fill Gaps · Mark Background Compounds 5. Peak Area Refinement - Compound Identification · Pathway Mapping · Fill Gaps · Mark Background Compounds 5. Peak Area Refinement - NoneMark Background Compounds Node- Unknown CompoundsSee "Marking Background Compounds in an Untargeted Analysis" on page 28.Normalize Areas NodeFill Gaps 6. Compound Identification - None 7. Compound Annotations Node · Group CompoundsNoneSearch Chemspider NodeGroup Compounds · Group Expected Compounds · Group Expected CompoundsSearch Mass Lists Node· Group Expected Compounds · Group CompoundsSearch mzCloud Node · Group CompoundsNoneSearch mzVault Node· Group Expected Compounds · Group Compounds 7. Pathway Mapping - Search Mass Lists · Search Mass Lists · Search ChemSpider · Map to KEGG Pathways · Map to BioCyc Pathways ·	Workflow node	Input nodes	Output nodes
Initial CompoundsMark Background CompoundsGroup Compounds NodeDetect Compounds· Compound Identification · Pathway Mapping · Fill Gaps5. Peak Area Refinement	Detect Compounds Node	Spectrum Processing	
Image: Parkway Mapping • Fill GapsParkway Mapping • Fill Gaps5. Peak Area RefinementNoneMark Background Compounds Node- Unknown CompoundsSee "Marking Background Compounds in an Untargeted Analysis" on page 28.NoneNormalize Areas NodeFill GapsMark Background Compounds6. Compound IdentificationKer Parking CompoundsNone7. Pedict Compositions NodeGroup CompoundsNone9. Predict Compositions NodeGroup CompoundsNone9. Search Mass Lists Node• Group Expected CompoundsNone9. Search Mass Lists Node• Group Expected CompoundsNone9. Search mzCloud Node• Group Expected CompoundsNone9. Search mzVault Node• Group CompoundsNone9. T. Pathway MappingAllGroup CompoundsNone9. Rompound Scoring· Search Mass Lists · Search Mass Lists · Map to KEGG Pathways · Map to Metabolika Pathways 	Fill Gaps Node	Group Compounds	
NoneSee "Marking Background Compounds in an Untargeted Analysis" on page 28.NoneNormalize Areas NodeFill GapsMark Background Compounds6. Compound IdentificationAssign Compound Annotations NodeGroup CompoundsNonePredict Compositions NodeGroup CompoundsNoneSearch ChemSpider NodeGroup CompoundsNoneSearch Mass Lists Node• Group CompoundsNoneSearch Mass Lists Node• Group Expected CompoundsNoneSearch mzCloud Node• Group Expected CompoundsNoneSearch mzVault Node• Group Expected CompoundsNoneSearch mzVault Node• Group CompoundsNoneAllGroup CompoundsNone8. Compound ScoringNoneApply Spectral Distance Node• Search Mass Lists • Map to KEGG Pathways • Map to BioCyc PathwaysNone	Group Compounds Node	Detect Compounds	Pathway MappingFill Gaps
Unknown CompoundsCompounds in an Untargeted Analysis" on page 28.Normalize Areas NodeFill GapsMark Background Compounds6. Compound IdentificationSearch ChemSpider NodeGroup CompoundsNonePredict Compositions NodeGroup CompoundsNoneSearch ChemSpider NodeGroup CompoundsNoneSearch Mass Lists Node• Group Expected CompoundsNoneSearch mzCloud Node• Group Expected CompoundsNoneSearch mzCloud Node• Group Expected CompoundsNoneSearch mzVault Node• Group Expected CompoundsNoneSearch mzVault Node• Group CompoundsNoneSearch mzVault Node• Group CompoundsNoneAllGroup CompoundsNone8. Compound Secring• Search Mass Lists • Search ChemSpider • Map to KEGG Pathways • Map to BioCyc Pathways • Map to BioCyc Pathways • Map to BioCyc Pathways • Map to BioCyc PathwaysNone	5. Peak Area Refinement		
6. Compound IdentificationI0IAssign Compound Annotations NodeGroup CompoundsNoneIPredict Compositions NodeGroup CompoundsNoneISearch ChemSpider NodeGroup CompoundsNoneISearch Mass Lists Node• Group Expected Compounds • Group CompoundsNoneISearch mzCloud Node• Group Expected Compounds • Group CompoundsNoneISearch mzVault Node• Group Expected Compounds • Group CompoundsNoneISearch mzVault Node• Group Compounds • Group CompoundsNoneIAllGroup CompoundsNoneI8. Compound Scoring• Search Mass Lists • Search ChemSpider • Map to KEGG Pathways • Map to BioCyc PathwaysNoneApply mzLogic Node• Search Mass Lists • Map to BioCyc Pathways • Map to BioCyc Pathways • Map to BioCyc Pathways • Map to BioCyc Pathways • Map to BioCyc PathwaysNone		Compounds in an Untargeted	None
Assign Compound Annotations NodeGroup CompoundsNonePredict Compositions NodeGroup CompoundsNoneSearch ChemSpider NodeGroup CompoundsNoneSearch Mass Lists Node- Group Expected CompoundsNoneSearch Mass Lists Node- Group Expected CompoundsNoneSearch mzCloud Node- Group Expected CompoundsNoneSearch mzVault Node- Group Expected CompoundsNoneSearch mzVault Node- Group CompoundsNoneJ. Pathway Mapping- Group CompoundsNoneAllGroup CompoundsNone8. Compound Scoring- Search Mass Lists - Search ChemSpider - Map to KEGG Pathways - Map to Metabolika Pathways - Map	Normalize Areas Node	Fill Gaps	Mark Background Compounds
Predict Compositions NodeGroup CompoundsNoneSearch ChemSpider NodeGroup CompoundsNoneSearch Mass Lists Node• Group Expected CompoundsNoneSearch Mass Lists Node• Group Expected CompoundsNoneSearch mzCloud Node• Group Expected CompoundsNoneSearch mzVault Node• Group Expected CompoundsNoneSearch mzVault Node• Group Expected CompoundsNoneT. Pathway MappingImage: CompoundsNoneAllGroup CompoundsNone8. Compound ScoringImage: CompoundsNoneApply Spectral Distance Node• Search Mass Lists · Search ChemSpider · Map to KEGG Pathways · Map to BioCyc PathwaysNoneApply mzLogic Node• Search Mass Lists · Search ChemSpider · Map to Metabolika Pathways · Map to Metabolika Pathways · Map to BioCyc PathwaysNone	6. Compound Identification		
Search ChemSpider NodeGroup CompoundsNoneSearch Mass Lists Node• Group Expected CompoundsNoneSearch Mass Lists Node• Group Expected CompoundsNoneSearch mzCloud Node• Group Expected CompoundsNoneSearch mzVault Node• Group Expected CompoundsNoneSearch mzVault Node• Group Expected CompoundsNoneSearch mzVault Node• Group CompoundsNoneSearch mzVault Node• Group CompoundsNoneAllGroup CompoundsNone8. Compound ScoringNoneApply Spectral Distance Node• Search Mass Lists • Search ChemSpider • Map to KEGG Pathways • Map to BioCyc Pathways • Map to BioCyc Pathways • Map to Metabolika Pathways • Map to Metabolika Pathways • Map to Metabolika Pathways • Map to Metabolika Pathways • Map to BioCyc Pathways • Map to BioCyc Pathways • Map to BioCyc PathwaysNone	Assign Compound Annotations Node	Group Compounds	None
Search Mass Lists NodeGroup Expected CompoundsNoneSearch mzCloud NodeGroup Expected CompoundsNoneSearch mzVault NodeGroup CompoundsNoneSearch mzVault NodeGroup Expected CompoundsNoneJ. Pathway MappingGroup CompoundsNoneAllGroup CompoundsNone8. Compound ScoringSearch Mass ListsNoneApply Spectral Distance NodeSearch Mass ListsNoneSearch ChemSpiderMap to KEGG PathwaysNoneApply mzLogic NodeSearch Mass ListsNoneMap to Metabolika PathwaysMap to Metabolika PathwaysNoneMap to BioCyc PathwaysMap to Metabolika PathwaysNoneMap to BioCyc PathwaysMap to Metabolika PathwaysNoneSearch ChemSpiderMap to Metabolika PathwaysNoneApply mzLogic NodeSearch Mass ListsNoneMap to BioCyc PathwaysMap to BioCyc PathwaysNoneSearch ChemSpiderMap to BioCyc PathwaysNoneSearch ChemSpiderMap to BioCyc PathwaysNone	Predict Compositions Node	Group Compounds	None
Search mzCloud NodeGroup Expected CompoundsNoneSearch mzVault NodeGroup CompoundsNoneSearch mzVault NodeGroup Expected CompoundsNone7. Pathway MappingGroup CompoundsNoneAllGroup CompoundsNone8. Compound ScoringSearch Mass ListsNoneApply Spectral Distance NodeSearch Mass ListsNoneSearch ChemSpiderMap to KEGG PathwaysNoneApply mzLogic NodeSearch Mass ListsNoneApply mzLogic NodeSearch Mass ListsNoneMap to BioCyc PathwaysMap to BioCyc PathwaysNoneMap to BioCyc PathwaysMap to BioCyc PathwaysNone	Search ChemSpider Node	Group Compounds	None
Group CompoundsSearch mzVault NodeGroup Expected Compounds Group Compounds7. Pathway MappingAllGroup Compounds8. Compound ScoringApply Spectral Distance NodeSearch Mass Lists Search ChemSpider Map to KEGG Pathways Map to BioCyc Pathways Search ChemSpiderApply mzLogic NodeSearch Mass Lists Map to Metabolika Pathways Map to BioCyc Pathways Map to BioCyc Pathways	Search Mass Lists Node		None
I Group Compounds7. Pathway MappingAllGroup CompoundsNone8. Compound ScoringNoneApply Spectral Distance Node• Search Mass Lists • Search ChemSpider • Map to KEGG Pathways • Map to BioCyc PathwaysNoneApply mzLogic Node• Search Mass Lists • Search ChemSpider • Map to Metabolika Pathways • Map to BioCyc PathwaysNone	Search mzCloud Node		None
AllGroup CompoundsNone8. Compound Scoring•Search Mass ListsNoneApply Spectral Distance Node•Search Mass ListsNone•Search ChemSpider•Map to KEGG PathwaysNone•Map to Metabolika Pathways•Map to BioCyc PathwaysNoneApply mzLogic Node•Search Mass ListsNone•Search ChemSpider•None•Map to Metabolika PathwaysNone•Map to Metabolika PathwaysNone•Map to Metabolika PathwaysNone•Map to BioCyc PathwaysNone	Search mzVault Node		None
8. Compound Scoring Apply Spectral Distance Node • Search Mass Lists None • Search ChemSpider • Map to KEGG Pathways None • Map to Metabolika Pathways • Map to BioCyc Pathways None Apply mzLogic Node • Search Mass Lists None • Map to Metabolika Pathways • Map to BioCyc Pathways None • Map to Metabolika Pathways • Map to Metabolika Pathways None • Map to Metabolika Pathways • Map to Metabolika Pathways None	7. Pathway Mapping		
Apply Spectral Distance Node• Search Mass ListsNone• Search ChemSpider • Map to KEGG Pathways • Map to Metabolika Pathways • Map to BioCyc Pathways• NoneApply mzLogic Node• Search Mass Lists • Search ChemSpider • Map to Metabolika Pathways • Map to Metabolika Pathways • Search ChemSpider • Map to Metabolika Pathways • Map to Metabolika Pathways • Map to BioCyc Pathways	All	Group Compounds	None
 Search ChemSpider Map to KEGG Pathways Map to Metabolika Pathways Map to BioCyc Pathways Apply mzLogic Node Search Mass Lists Search ChemSpider Map to Metabolika Pathways Map to Metabolika Pathways Map to Metabolika Pathways Map to BioCyc Pathways 	8. Compound Scoring		
 Search ChemSpider Map to Metabolika Pathways Map to BioCyc Pathways 	Apply Spectral Distance Node	Search ChemSpiderMap to KEGG PathwaysMap to Metabolika Pathways	None
Calculate Mass Defect Node Group Compounds None	Apply mzLogic Node	Search ChemSpiderMap to Metabolika Pathways	None
	Calculate Mass Defect Node	Group Compounds	None

Table 28. Input and output nodes for each workflow node (Sheet 2 of 3)

Workflow node	Input nodes	Output nodes
Compound Class Scoring Node	Group Compounds	None
Generate Molecular Networks Node (Beta)	Assign Compound Annotations	None
Pattern Scoring Node	Group Compounds	None
9. Expected Compounds		
FISh Scoring Node	Group Expected Compounds	None
Find Expected Compounds Node	Generate Expected CompoundsSpectrum Processing	Group Expected CompoundsMerge Features
Generate Expected Compounds Node	None	Find Expected CompoundsFilter By Mass Defect
Group Expected Compounds Node	Find Expected Compounds	FISh ScoringMark Background Compounds
Mark Background Compounds Node– Expected Compounds	Group Expected Compounds	None
Merge Features Node	Find Expected CompoundsDetect Compounds	None
10. Post-Processing (white nodes)		
Descriptive Statistics Node	None	None
Differential Analysis Node	None	None
Export Xcalibur Inclusion or Exclusion List Node	None	None

Table 28. Input and output nodes for each workflow node (Sheet 3 of 3)

Saving a Processing Workflow as a Template

You can run an analysis without saving the processing workflow in the Workflow Tree pane; however, you might want to save the processing workflow to a template for reuse later on. When you save the processing workflow as a template, the application does not automatically store it in the study folder. You can save a processing template to the Common Templates folder or a folder of your choice.

Note When you run an analysis, the application automatically saves the processing workflow to the result file. Selecting a result file on the Analysis Results page of a study and clicking **Reprocess** opens the processing workflow saved with the result file.

* To save a processing workflow as a template

- 1. Do one of the following:
 - To save the template in the Common Templates folder, click Save Common.

The Save Workflow dialog box opens to the Common Templates folder.

• To save the template to the last folder that you opened or to another folder, click **Save**.

The Save Workflow dialog box opens to the last opened folder.

2. Select the folder where you want to store the template, name the template, and click Save.

If the processing workflow is valid, the application saves the template with the file name extension .cdProcessingWF. If the processing workflow contains an error, an error message box opens.

3. If the Exporting Template Workflow Failed message box opens, read the list of errors, close the message box, fix the errors, and click **Save**.

Use the Workflows page (in any study) to create or edit processing workflows. The Workflows page is a tabbed page to the right of the Grouping & Ratios page and is available only when the Analysis pane is open.

6 Creating and Editing Processing Workflows Saving a Processing Workflow as a Template

7

Workflow Nodes

A processing workflow consists of a set of connected workflow nodes and each workflow node has a set of associated parameters. To create and edit processing workflows, see Chapter 6, "Creating and Editing Processing Workflows."

Note The following nodes are new in Compound Discoverer 3.1—Apply Spectral Distance, Calculate Mass Defect, Generate Molecular Networks (Beta), and Scripting Node.

On the Workflows page, in the Workflow Nodes pane, the workflow nodes are organized into functional groups. For information about the workflow nodes, see the following topics.

Contents

- Input and Output Nodes
- Spectrum Processing Nodes
- Trace Creation Nodes
- Compound Detection Nodes
- Peak Area Refinement Nodes
- Compound Identification Nodes
- Pathway Mapping Nodes
- Compound Scoring Nodes
- Expected Compounds Nodes
- Post-Processing Nodes

Input and Output Nodes

These topics describe the input and output nodes:

- Export Spectra Node
- Input Files Node

Export Spectra Node

Use the Export Spectra node to export all or a subset of the mass spectrum scans in an Xcalibur RAW file to an open-source format file. The Export Spectra node does not export the data from analog detectors.

Note You can add multiple Export Spectra nodes to a processing workflow.

The Export Spectra node requires input from one of the data processing nodes. The processing workflow shown in Figure 69 reads the input files, extracts the MS scans of interest, and exports the data to an open source format file. Running this processing workflow does not require an active software license.

Figure 69. Minimum processing workflow for the Export Spectra node

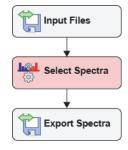


Table 29 describes the parameters for the Export Spectra node.

Parameter	Description
1. Output Data	
File Name	Specifies the file name of the exported file. If you leave this box empty, the node uses the result file name.
File Name Suffix	Specifies the suffix that the application appends to the file name of

the exported file.

Table 29. Export Spectra node parameters (Sheet 1 of 2)

Thermo Scientific

Parameter	Description
Export Format	Specifies the data format of the exported file.
	Selections:
	 DTA archive (*.dta.zip)—Generates a zip file that contains an individual DTA file for each MS scan in the raw data file. The individual file names include the scan number of the scan. The DTA file format is a simple text file that contains a peak table of mass and intensity. This format does not store information about data acquisition, such as the instrument method or mass resolution, and you cannot reprocess DTA files with a mass spectrometry application.
	• Mascot Generic Format (*.mgf)—Generates an MGF file, which lists the MS scans by retention time. The scan data for each time point consists of two columns: mass and intensity.
	• mzDATA (*mzData)—Generates an XML-based file that third-party mass spectrometry software packages can read.
	• mzML (*mzML)—Generates an XML-based file that third-party mass spectrometry software packages can read.

Table 29. Export Spectra node parameters (Sheet 2 of 2)

Input Files Node

Every processing workflow must begin with the Input Files node. This node has no parameters.

To view information about a result file's input files, open its Input Files Table.

Spectrum Processing Nodes

These nodes extract the mass spectral data from the input file set:

- Align Retention Times Node
- Filter By Mass Defect Node
- Filter By Scan Event Node
- Filter Centroids Node
- Select Spectra Node

Align Retention Times Node

Use the Align Retention Times node for chromatographic alignment of multiple input files. The Align Retention Times node compensates for small differences in the retention times of the components in a sequence of sample runs.

IMPORTANT In the spectrum processing node that supplies spectra to the Alignment, do not apply a scan filter that excessively reduces the retention time window, as doing so might cause an alignment failure. The Align Retention Times node requires a minimum amount of representative scan data to chromatographically align the input files in an analysis.

Table 30 describes the parameters for the Align Retention Times node.

Parameter	Description
1. General Settings	
Alignment Model	Specifies the curve fitting algorithm that the workflow uses to chromatographically align the input files.
	Default: Adaptive Curve
	Selections: Adaptive Curve or Linear
Alignment Fallback	Specifies the alternate model to apply when the Adaptive Curve regression model fails.
	Default: Use Linear Model
	Selections:
	• None—There is no alternative. Continue to use the Adaptive Curve model.
	 Don't Align—Do not chromatographically align the input files.
	• Use Linear Model—Use the Linear model instead of the Adaptive Curve model.
Maximum Shift [min]	Specifies the maximum retention time shift between the alignment features (chromatographic peaks with the same $m/z \times RT$ dimensions) in the input files.
	Default: 2 (± 2 minutes for each feature); range: 0.01–4.0

Table 30. Align Retention Times node parameters (Sheet 1 of 2)

Parameter	Description
Shift Reference File	Specifies whether to shift the retention time of all the detected features to eliminate any negative retention time values in the input file set.
	Default: True
	 Selections: True: If the data set includes features with negative retention times, the algorithm shifts all of the features to avoid cropping. False: Removes features with negative retention time values from the analysis (crops the feature from the feature list passed to the connected nodes).
Mass Tolerance	Specifies the mass tolerance to be used for feature matching.
	Default: 5.0 ppm; range: 0.1–50 ppm
Remove Outlier	Specifies whether the retention time algorithm ignores outlier landmark features.
	Default: True

Table 30. Align Retention Times node parameters (Sheet 2 of 2)

Related Topics

- Viewing the Corrected Retention Times of the Alignment Features
- File Alignments Table

Filter By Mass Defect Node

Use the Filter By Mass Defect node to keep or remove mass spectral peaks (centroids) in the full (MS1) scan data that fall within a set of specified mass tolerance and mass defect windows.

You can add multiple Filter By Mass Defect nodes to a processing workflow.

To specify the elemental compositions for the node, you can enter the elemental compositions in the Custom Compositions area of the node, or you can use the Generate Expected Compounds Node to provide the compositions. Table 31 describes the parameters for the Filter By Mass Defect node.

 Table 31.
 Filter By Mass Defect node parameters (Sheet 1 of 2)

Parameter	Description
1. General Settings	
Filter Direction	Specifies whether the selected mass defect filter keeps or removes mass spectral peaks (centroids) from further processing.
	Default: Keep Selections: Keep or Remove
Mass Defect Type	Specifies the mass defect type.
	Default: Standard Mass Defect
	Selections:
	Fractional Mass = exact mass – floor (exact mass)
	Standard Mass Defect = exact mass – nominal mass
	Relative Mass Defect = $\frac{1e6 \times (exact mass - nominal mass)}{(exact mass)}$
	Kendrick Mass Defect = Kendrick mass – nominal Kendrick mass
	Where:
	Exact mass = Monoisotopic mass of the elemental composition
	Nominal mass = Integer mass
	Calculates the integer mass using the selected rounding function (floor, ceiling, or round)
	Kendrick Mass = exact mass × nominal mass of Kendrick formula exact mass of Kendrick formula
Kendrick Formula	When you select the Kendrick Mass Defect type, this user-specified elemental composition specifies the Kendrick formula.
Nominal Mass Rounding	Specifies how the node calculates nominal masses.
	Default: Floor
	Selections:
	• Floor rounds down.
	• Ceiling rounds up.
	 Round rounds to the nearest integer value.

i arameter	Description
2. Tolerances	
	fect values calculated above and the exact mass values calculated from the ition input, these mass tolerance values define the rectangular mass defect
Mass Tolerance	The input from the data processing nodes is a table of m/z values and intensities for each full (MS1) scan.
	Specifies the mass tolerance for the ions that the filter removes or passes through to the next node.
	Default: 50 Da
	Selection: 0 to 6000 Da
Mass Defect Tolerance	Specifies the mass defect tolerance for the ions that the filter removes or passes through to the next node.
	Default: 0.025 (Da or unit-less for the Relative Mass Defect selection)
	Range: 0–No limit
3. Custom Composi	tions
Composition (5 entry boxes)	Specify the elemental compositions that the node uses to create the mass defect filters.
	Leave these boxes empty if you want to use one or more Generate Expected Compounds nodes to generate the list of elemental compositions.
Ions	Specifies the ion definitions to be used with the custom compositions. For each elemental composition, the node creates one mass defect filter for each ion definition—that is, if you select five ion definitions, the node creates five mass defect filters for each elemental composition.

Select the ion definitions from the dropdown list. Use the Ion

Definition Editor to create new ion definitions.

Table 31. Filter By Mass Defect node parameters (Sheet 2 of 2)

Description

Related Topics

Parameter

- Adding or Editing Ion Definitions with the Ion Definition Editor
- Calculating the Mass Defect of an Elemental Composition

Filter By Scan Event Node

Use the Filter By Scan Event node to filter the mass spectra by scan events.

Table 32 describes the parameters for the Filter By Scan Event node.

 Table 32.
 Filter By Scan Event node parameters (Sheet 1 of 3)

Parameter	Description
Filter Settings	
Mass Analyzer	Specifies the mass analyzer that the instrument used to acquire the scan. A Thermo Scientific hybrid mass spectrometer, such as the LTQ Orbitrap mass spectrometer, contains two mass analyzers and can acquire both ITMS (low-resolution) and FTMS (high-resolution) scans in one data file.
	Filter selection: Any, Is, Is Not Default: (Not Specified)—The application does not filter scan events by the mass analyzer that was used to acquire the data.
	Check box selections: • Ion Trap (ITMS) • Fourier Transform (FTMS)) • Time of Flight (TOFMS) • Single Quad (SQMS) • Triple Quad (TSMS) • Sector Field (SectorMS)
MS Order	Specifies the MS order (scan power that the instrument used) of the scans that you want the node to filter. Filter selection: Any, Is, Is Not Default: (Not Specified)—The node does not filter the scans by MS order.
	Check box selections: MS1–MS7
	Note The Detect Compounds and Find Expected Compounds nodes search the full (MS1) scans for mass peaks. If you filter out the MS1 scans by selecting Is MS2 or higher for the MS Order, the result tables for these nodes are empty.

Parameter	Description
Activation Type	Specifies the activation types that the instrument used to produce the product ion spectra.
	Filter selection: Any, Is, Is Not Default: (Not Specified)—The node does not filter the scans by activation type.
	 Check box selections: CID (Collision Induced Dissociation) MPD (Multi Photon Dissociation) ECD (Electron Capture Dissociation) PQD (Pulsed Q Collision Induced Dissociation) ETD (Electron Transfer Dissociation) HCD (Higher Energy Collision Dissociation) EThcD (ETD With Supplemental HCD) UVPD (Ultra Violet Photo Dissociation) Any Activation Type
Min. Collision Energy	Specifies the minimum Normalized Collision Energy [™] for a higher-order scan to pass through the filter. Default: 0 (no filtering) Minimum value: 0 Maximum value: Unchecked
Max. Collision Energy	Specifies the maximum Normalized Collision Energy for a higher-order scan to pass through the filter. Default: 1000 Minimum value: 0 Maximum value: 1000

 Table 32.
 Filter By Scan Event node parameters (Sheet 2 of 3)

Parameter	Description
Scan Type	Specifies the scan type for the scan event that the instrument used to produce the product ion.
	Filter selection: Any, Is, Is Not Default: (Not Specified)—The node does not filter the scan event by scan type.
	Check box selections: • Full • Single Ion Monitoring (SIM) • Single Reaction Monitoring (SRM)
Polarity Mode	Specifies the polarity mode for the scan.
	Filter selection: Any, Is, Is Not
	Default: (Not Specified)—The node does not filter the scan event by polarity mode.
	Check box selections: • Positive • Negative

 Table 32.
 Filter By Scan Event node parameters (Sheet 3 of 3)

Filter Centroids Node

Use the Filter Centroids node to remove mass spectral peaks (centroids) that are below a user-specified intensity threshold from the mass scans, below a user-specified signal-to-noise threshold for FTMS scans, or both.

Table 33 describes the parameters for the Filter Centroids node.

Parameter	Description
1. General	
S/N Threshold (for FT-only)	Specifies the minimum signal-to-noise threshold for each centroid in an FTMS scan. The node excludes centroids from the analysis that are below this intensity value.
	The application uses the spectrum noise reported by an instrument.
	Default: 1.5
Minimum Intensity Threshold	Specifies the minimum intensity threshold for the mass spectral peaks (centroids). The node excludes centroids from the analysis that are below this intensity value.
	Default: 0 (no filtering)

Select Spectra Node

The raw data file (Xcalibur RAW file) contains the mass spectral scans acquired by your Thermo Scientific mass spectrometer and any optional data acquired by a PDA, UV-VIS, or analog detector during the acquisition run. The Select Spectra node can read and filter the mass spectral scan data. The Select Spectra node cannot read the optional data acquired by a PDA, UV-VIS, or analog detector.

The mass spectral scans are numbered 1, 2, 3, and so on from the beginning to the end of the acquisition run. Use the Select Spectra node to select the scans that you want the application to process. Limiting processing to the scans of interest decreases processing time and minimizes false positives. For example, if you know the retention time of the compounds of interest, exclude scans that fall outside a specific retention time window.

IMPORTANT When using settings other than the defaults for the Select Spectra node, follow these guidelines:

- The application uses the full (MS1) scans to measure the accurate mass and isotope patterns of the mass spectral peaks; therefore, do not filter out the full (MS1) scans when the processing workflow includes the Find Expected Compounds and Detect Compounds nodes.
- Both the Align Retention Times and Find Expected Compounds nodes require a representative amount of data to function properly. If you excessively reduce the retention time window (for example, by using an RT or scan number range), the Align Retention Times node's chromatographic alignment algorithm and the Find Expected Compounds node's automatic peak width detection algorithm might fail to produce satisfactory results.

Table 34 describes the parameters in the Select Spectra node.

Parameter	Description
1. General Settings	
Precursor Selection	Specifies the MS order of the precursor scans for higher-order MS ⁿ scans, like MS ³ , MS ⁴ , and so forth.
	Default: Use MS(n – 1) Precursor
	Selections: Use MS1 Precursor or Use MS(n - 1) Precursor

Table 34. Select Spectra node parameters (Sheet 1 of 8)

Parameter	Description
Use Isotope Pattern in Precursor Reevaluation	Determines whether the node considers the isotope pattern in reevaluating precursors.
	Default: True Selections:
	 True—The node considers the isotope pattern in reevaluating precursors.
	• False—The node does not consider the isotope pattern in reevaluating precursors.
Provide Profile Spectra	When set to True, the node stores the profile data for the scans.
	When set to Automatic, the node checks whether any other nodes in the processing workflow that require profile data. If the profile data is not needed, the node only stores the centroid data.
	Default: Automatic
Store Chromatograms	Specifies whether the application creates a TIC trace and a BPC trace for each input file and lists these traces in the Specialized Traces table.
	Default: False

Table 34. Select Spectra node parameters (Sheet 2 of 8)

2. Spectrum Properties Filter

Note The retention time filter excludes scans outside the specified limits; however, the application does not check the validity of the retention time settings against the actual acquisition time for the raw data file. When both the lower and upper RT limits are set to 0 (default), the application does not use retention time to filter scans.

Lower RT Limit	Excludes scans acquired before the user-specified retention time.
	Default: 0
	Minimum value: 0
	Maximum value: Unchecked
Upper RT Limit	With the exception of a setting of 0, excludes scans that were acquired after the user-specified retention time.
	Default: 0; Minimum value: 0; Maximum value: Unchecked
	ber filter excludes scans outside the specified limits; however, the

application does not check the validity of the scan number settings against the actual scan numbers in the raw data file. When both the first and last scan number are set to 0 (default), the application does not filter scans by scan number. When filtering by scan number, verify the scan number range in a mass spectrometry viewer application, such as Qual Browser or FreeStyle.

Parameter	Description
First Scan	Specifies the scan number of the first available scan that you wan the node to process.
	When this parameter is set to 0, the node processes the first scar that passes through the other filters.
	When this parameter is set to a value that is greater than the last available scan number, the node filters out all of the scans.
	Default: 0 (no filtering)
	Minimum value: 0; Maximum value: Unchecked
Last Scan	Specifies the last available scan number that you want the node process.
	When this parameter is set to 0 or a value that is greater than the last available scan number, the node processes the last scan that passes through the other filters.
	Default: 0 (no filtering)
	Minimum value: 0; Maximum value: Unchecked
Ignore Specified Scans	Specifies the scan numbers that the analysis ignores. Placing the cursor in the Ignore Specified Scans box and clicking the More icon,, opens the Edit Parameter Text for Ignore Specified Scans dialog box where you can do the following:
	• Manually enter the scan numbers to ignore.
	• Load a text file that lists the scans numbers to ignore.
	• Create and save a list of scans numbers to ignore.
	Edit parameter text for Ignore Specified Scans
	Enter single scan numbers or scan ranges (e.g., 123-234, or 123-, or - +123) on separate lines.

Table 34. Select Spectra node parameters (Sheet 3 of 8)

Parameter	Description
U	filter excludes higher-order scans of precursor ions with a charge specified limits. The charge state filter does not affect the MS1
Lowest Charge State	Excludes higher-order scans of precursor ions with a lower charge state than the specified charge state.
	Default: 0; Minimum: 0
Highest Charge State	Filters out scans from precursor ions with a higher charge state than the specified charge state.
	Default: 0 (specifies no upper limit)
Min. Precursor Mass	Specifies the minimum precursor mass for a higher-order scan.
	Default: 0 (no filtering)
	Minimum value: 0; Maximum value: Unchecked
Max. Precursor Mass	Specifies the maximum precursor mass for a higher-order scan.
	Default: 0 (no filtering)
	Minimum value: 0; Maximum value: Unchecked
Total Intensity Threshold	Excludes scans that fall below the specified total intensity threshold. The total intensity of a mass spectrum is the summed intensity of its mass spectrum peaks (centroids).
Minimum Peak Count	Specifies the minimum number of mass spectrum peaks (centroids) that must be in the spectrum for the scan to pass through the filter.
	Minimum value: 1; Maximum value: Unchecked

 Table 34.
 Select Spectra node parameters (Sheet 4 of 8)

Parameter	Description
3. Scan Event Filters	
Mass Analyzer	Specifies the mass analyzer that the instrument used to acquire the scan. An LTQ Orbitrap hybrid mass spectrometer can acquire both ITMS (low-resolution) and FTMS (high-resolution) scans in one data file.
	Filter selection: Any, Is, Is Not Default: (Not Specified)—The node does not filter scan events by the mass analyzer used to acquire the data.
	Check box selections: Ion Trap (ITMS) Fourier Transform (FTMS)) Time of Flight (TOFMS) Single Quad (SQMS) Triple Quad (TSMS) Sector Field (SectorMS)
MS Order	Specifies the MS order of the scans that you want the node to process.
	Filter selection: Any, Is, Is Not Default: Any—The application does not filter the scans by MS order.
	Check box selections: MS1–MS7
	Note The Detect Compounds and Find Expected Compounds nodes search the MS1 scans for mass peaks. If you filter out the MS1 scans by selecting Is MS2 or higher for the MS Order, the result tables for these nodes are empty.

Table 34.Select Spectra node parameters (Sheet 5 of 8)

Parameter	Description
Activation Type	Specifies the activation type that the instrument used to produce the product ion spectra.
	Filter selection: Any, Is, Is Not Default: (Not Specified)—The node does not filter the scans by activation type.
	 Check box selections: CID (Collision Induced Dissociation) MPD (Multi Photon Dissociation) ECD (Electron Capture Dissociation) PQD (Pulsed Q Collision Induced Dissociation) ETD (Electron Transfer Dissociation) HCD (Higher Energy Collision Dissociation) EThcD (ETD With Supplemental HCD) UVPD (Ultra Violet Photon Dissociation) Any Activation Type
Min. Collision Energy	Specifies the minimum Normalized Collision Energy for a higher-order scan to pass through the filter. Default: 0 (no filtering)
	Minimum value: 0; Maximum value: Unchecked
Max. Collision Energy	Specifies the maximum Normalized Collision Energy for a higher-order scan to pass through the filter.
	Default: 1000; Minimum value: 0; Maximum value: 1000
Scan Type	Specifies the scan type for the scan event that the instrument used to produce the product ion.
	Filter selection: Any, Is, Is Not Default: Any—The application does not filter the scan events by scan type.
	 Check box selections: Full Selected Ion Monitoring (SIM) Selected Reaction Monitoring (SRM)

 Table 34.
 Select Spectra node parameters (Sheet 6 of 8)

Parameter	Description
Polarity Mode	Specifies the polarity mode for the scan.
	Filter selection: Any, Is, Is Not
	Default: (Not Specified)—The node does not filter the scan events by polarity mode.
	Check box selections: • Positive • Negative
4. Peak Filters	
S/N Threshold (FT-only)	Specifies the signal-to-noise threshold for mass peaks in an FTMS scan. Mass peaks below this threshold are filtered out.
	Default: 1.5
5. Replacements for Unre	cognized Properties
Unrecognized Charge Replacements	Specifies the charge state or states to process when the charge state of the precursor ion is indeterminate.
	Default: 1
	In the Qual Browser window, an indeterminate charge state is specified with a question mark label (z=?) in a spectrum cell. In the FreeStyle application, ions with indeterminate charge states are labeled with a charge state of 0.
	Default: 1; Selections: All, 1, 2, 3, 4, 5, 6, 7, and 8
Unrecognized Mass Analyzer Replacements	Specifies the mass analyzer used to acquire the data when the application cannot retrieve this information from the input file.
	Default: ITMS
	Selections: Ion Trap (ITMS) Fourier Transform (FTMS) Time of Flight (TOFMS) Single Quad (SQMS) Triple Quad (TSMS) Sector Field (SectorMS)
Unrecognized MS Order Replacements	Specifies the MS order when the application cannot retrieve this information from the input file.
	Default: MS2; Selections: MS1–MS10

Table 34. Select Spectra node parameters (Sheet 7 of 8)

Parameter	Description
Unrecognized Activation Type Replacements	Specifies the activation type when the application cannot retrieve this information from the input file.
	Default: CID
	 Selections: CID (Collision Induced Dissociation) MPD (Multi Photon Dissociation) ECD (Electron Capture Dissociation) PQD (Pulsed Q Collision Induced Dissociation) ETD (Electron Transfer Dissociation) HCD (High Energy Collision Dissociation) EThcD (ETD With Supplemental HCD) UVPD (Ultra Violet Photon Dissociation)
Unrecognized Polarity Replacement	Specifies the polarity mode when the application cannot retrieve this information from the input file. Default: (+) Selections: Positive (+) or Negative (-)
Unrecognized MS Resolution @ 200 Replacements	Specifies the resolution at m/z 200 for MS scans when the node cannot retrieve the resolution from the scan header.
	Default: 60 000
Unrecognized MSn Resolution @ 200	Specifies the resolution at m/z 200 for MS/MS scans when the node cannot retrieve the resolution from the scan header.
Replacements	Default: 30 000

Table 34. Select Spectra node parameters (Sheet 8 of 8)

Trace Creation Nodes

These nodes create specialized chromatographic traces:

- Create Analog Trace Node
- Create FISh Trace Node
- Create Mass Trace Node
- Create Pattern Trace Node

Create Analog Trace Node

Use the Create Analog Trace node to view the chromatograms for these trace types: ultraviolet-visible (UV), photo-diode array (PDA), or analog. Your Thermo Scientific data system supports several brands of UV-Vis and PDA detectors. You can acquire UV traces from a UV-Vis or a PDA detector and PDA traces from a PDA detector. If your analog detector is not supported by a Thermo Scientific data system, you can acquire an analog trace by connecting the detector to one of the analog channels on the communications panel of your Thermo Scientific mass spectrometer.

You can access analog traces from the Specialized Traces Table.

Note The Create Analog Trace node can also convert and display a pressure trace from an LC pump and a temperature trace from a column heater or autosampler with temperature control, when these instruments are controlled by the Xcalibur data system or equivalent Thermo Scientific application.

Table 35 describes the parameters for the Analog Traces node.

Parameter	Description
1. General Settings	
Trace Type	Specifies whether the application extracts a UV, PDA, or analog trace from the raw data file.
	Default: UV; Selections: UV, PDA, or Analog
RT Offset [min]	Specifies the offset time, in minutes, between the UV, PDA, or analog detector and the mass spectrometer traces.
	A negative value shortens and a positive value lengthens the apparent retention time of the peaks detected by the UV, PDA, or analog detector.
	Default: 0; range: –10 to 10
Custom Label	Type text to identify the trace in the Specialized Traces table on the result file page.
2. PDA Settings	
Total Scan	Specifies whether the node extracts a total scan trace for the scanned wavelength range.
	Total scan traces display the average absorbance for each time point of all the wavelengths in the scan range.
	Default: False

 Table 35.
 Create Analog Trace node parameters (Sheet 1 of 2)

Parameter	Description
Spectrum Maximum	Specifies whether the node extracts a spectrum maximum trace for the scanned wavelength range.
	Spectrum maximum traces display a plot of the maximum absorbance values in the scan range for each time point.
	Default: False
Note Use the Min. ar versus time.	nd Max. Wavelength boxes to specify a trace of average absorbance
	romatogram for a specific scan wavelength, type the same per in the Min. and Max. Wavelength boxes.
beginning waveler	of the average absorbance values for a range of wavelengths, type the ngth number in the Min. Wavelength box and the ending her in the Max. Wavelength box.
Wavelength Range	Specifies whether the node extracts the entire acquired scan or a wavelength range.
	Default: True—Uses a specified wavelength range.
Min. Wavelength	Default: Irue—Uses a specified wavelength range. Specifies the beginning wavelength, in nanometers, of the trace that you want the node to extract.
Min. Wavelength	Specifies the beginning wavelength, in nanometers, of the trace
Min. Wavelength Max. Wavelength	Specifies the beginning wavelength, in nanometers, of the trace that you want the node to extract.

Table 35. Create Analog Trace node parameters (Sheet 2 of 2)

Create FISh Trace Node

Use the Create (Fragment Ion Search) FISh Trace node to create FISh traces. The Create FISh Trace node requires fragmentation scans and takes input from any of the data processing nodes. The output from the FISh trace node is a summed FISh trace that is accessible from the Specialized Traces Table, individual fragment traces that are accessible from the FISh Trace Fragments result table, or both.

Table 36 describes the parameters for the Create FISh Trace node.

Parameter	Description	
1. Compound Selection		
Compound	Specifies the compound that the node uses to generate expected fragment ions.	
	The selection list contains the compounds in the user-created Expected Compounds library .	
	IMPORTANT To run an analysis that includes the Create FISh Trace node, you must select a compound from this list.	
2. Trace Settings		
Mass Tolerance	Specifies the mass tolerance that the node uses to create the FISh trace.	
	Default: 2.5 mmu	
Summed Trace	Specifies whether the node generates a summed trace of all the detected fragment ions. You can access the summed FISh trace from the Specialized Traces result table.	
	Default: True	
Individual Traces	Specifies whether the node generates individual traces for each generated fragment ion. You can access the individual traces from the FISh Trace Fragments result table.	
	Default: True	
Custom Label	Type text that you can use to identify the chromatogram in a report.	
	This box accepts alphanumeric and special characters.	
3. Scan Filter Settings		
Scan Polarity	Specifies the polarity of the scan.	
	Default: + (Positive); Selections: Positive or Negative	

Table 36. Create FISh Trace node parameters (Sheet 1 of 3)

Parameter	Description
Fragment Mode	Specifies the fragmentation mode of the fragmentation scans that you want the node to extract.
	Default: Data-Dependent; selections: Data-Dependent and Data Independent
	Select Data-Dependent for data-dependent fragmentation (DDF scans or Data Independent for all-ion fragmentation (AIF) scans
4. Fragment Prediction	Settings
Use General Rules	Specifies whether the node uses the general fragmentation rules fo fragment prediction.
Use Libraries	Specifies whether the node uses the fragmentation libraries for fragment prediction.
	Default: True
Max. Depth	Specifies the maximum number of steps in the fragmentation pathway.
	Range: 1 to 20
Aromatic Cleavage	Specifies whether the node includes a cleavage step in the fragmentation pathway for highly aromatic structures—that is, fo aromatic structures where <i>n</i> in Huckel's rule is 2 or higher.
	4n + 2 = 0 or a positive integer for the number of electrons in the delocalized, conjugated p-orbital cloud
	For example, the following structure is aromatic with an n value of 2.
	HON
	Default: True

 Table 36.
 Create FISh Trace node parameters (Sheet 2 of 3)

Parameter	Description
Min. Fragment <i>m</i> /z	Specifies the minimum m/z value of a fragment ion to be generated by the prediction fragmentation pathway.
	Default: 50; range: 0 or higher
Max. Fragment <i>m</i> / <i>z</i>	Specifies the maximum <i>m/z</i> value of a fragment ion to be generated by the prediction fragmentation pathway.
	When the value is set to 0, the node ignores this parameter.
	Default: 0

Table 36. Create FISh Trace node parameters (Sheet 3 of 3)

Related Topics

- Modifying the Expected Compounds List
- Troubleshooting Common Analysis Errors

Create Mass Trace Node

Use the Create Mass Trace node to extract a mass chromatogram that you can access from the Specialized Traces Table of the result file. You can specify the type, the fragmentation order, and the polarity of the trace. For an XIC trace, you must specify the mass range.

Table 37 describes the parameters for the Create Mass Trace node.

Table 37. Create Mass Trace node parameters (Sheet 1 of 2)

Parameter	Description
1. General Settings	
Trace Type	Specifies the chromatogram type to be generated.
	Default: BPC
	Selections: TIC (total ion chromatogram), BPC (base peak chromatogram), or XIC (extracted ion chromatogram)
MS Order	Specifies the MS order of the mass spectra that make up the chromatogram.
	Default: MS1; Selections: MS1–MS10
Polarity	Specifies the ionization polarity used to produce the mass spectra that make up the chromatogram.
	Default: + (Positive); Selections: + (Positive) or – (Negative)

Parameter	Description
Custom Label	Use this box to type text that you can use to identify the chromatogram.
	This text appears in the Custom Label column of the Specialized Traces result table.
	This box accepts alphanumeric and special characters.
2. XIC Settings	
Mass [Da]	Defines the mass-to-charge (m/z) value of the extracted ion chromatogram (XIC).
	Default: 0
Mass Tolerance	Specifies the mass tolerance for the spectral search.
(typed numeric value and selected units)	When you select Da or mmu (0.001 Da) in the units list, the mass tolerance is an absolute \pm value for the mass specified in the Mass box.
	When you select ppm (parts per million) in the Units list, the mass tolerance is a relative range:
	Mass ± (Mass × User-specified ppm)/1e6
	Default: 3 ppm
	Range: 0 to no upper limit; units: Da, mmu, ppm

Table 37. Create Mass Trace node parameters (Sheet 2 of 2)

Create Pattern Trace Node

Use the Create Pattern Trace node to draw a chromatogram from the mass peaks that match a specific pattern within the filtered set of spectra. The pattern can be based on the elemental composition of a target compound or on a user-specified pattern. To view the pattern trace in the Chromatogram view for a result file, open the Specialized Traces Table and select the pattern trace.

Table 38 describes the parameters for the Create Pattern Trace node.

Parameter	Description		
1. General Settings			
Isotope Ratios	Displays either the elemental composition of the compound of interest or the following text: [custom pattern].		
	You use the Isotope Ratio Editor to set up the isotope pattern for the compound of interest.		
	The application can automatically set up the isotope pattern for a non-isotopically labeled compound by using its elemental composition. For isotopically-labeled compounds, you must enter the expected mass shifts of the isotopic peaks as well as their relative intensity to the A0 isotope.		
	IMPORTANT In the defined processing workflows, this parameter is set to C15S with the following three isotope selections:		
	• Monoisotopic ion (100% intensity)		
	• A2 ion with one sulfur-34 atom (peak with a +1.9958 Da mass shift and a 4.52% relative intensity)		
	• A2 ion with two carbon-13 atoms (peak with a +2.00669 Da mass shift and a 1.27% relative intensity)		
	When you create a new processing workflow, you must specify the isotope ratios of interest to run an analysis.		
Mass Tolerance	Specifies the mass tolerance for the mass shifts between the mass spectral peaks in the pattern.		
	Range: 0.0–1e6 ppm; default: 5 ppm		
	You set up the pattern with the Isotope Ratio Editor.		
Intensity Tolerance [%]	Specifies the relative intensity tolerance of the mass spectral peaks in the pattern.		
	Range: 0.01–100.0		
	The A0 isotope is always the isotope with the lowest m/z value, but it is not necessarily the isotope with the highest intensity. For example, with more than one bromine atom, a bromine and a chlorine atom, or more than four chlorine atoms, the M + 2 (A2) isotope is the most intense isotope.		

 Table 38.
 Create Pattern Trace node parameters (Sheet 1 of 2)

Parameter	Description	
MS Order	Specifies the MS order of the mass spectrum.	
	Selections: MS1–MS10	
Polarity	Specifies the polarity of the mass spectrum.	
	Selections: Positive or Negative	
Custom Label	Use this box to enter a description of the trace.	

Table 38. Create Pattern Trace node parameters (Sheet 2 of 2)

Related Topics

- Setting Up Isotope Patterns for Pattern Scoring or Pattern Traces
- Specialized Traces Table

Compound Detection Nodes

Use these nodes to detect unknown compounds:

- Analyze Labeled Compounds Node
- Detect Compounds Node
- Fill Gaps Node
- Group Compounds Node

Analyze Labeled Compounds Node

Use the Analyze Labeled Compounds node to detect labeled compounds and their labeling rates.

Table 39.	Analyze L	abeled Com	pounds node	(Sheet 1	of 3)
-----------	-----------	------------	-------------	----------	-------

Parameter	Description	
1. Label Settings		
Label Element	Specifies the labeled element and its isotope in this format:	
	[mass number]element symbol	
	Where <i>element symbol</i> is the symbol in the periodic table	

, D	
Parameter	Description
Max. Exchange	Specifies the maximum number of exchangeable atoms for any of the compounds. If the number of exchangeable atoms for a compound is below this number, the application uses the lower number.
	If set to zero, the compound's elemental composition determines the maximum number of exchangeable atoms.
	Default: 25
Source Efficiency [%]	Specifies the isotopic purity of the labeled compound introduced into the biological system.
	Most commercially available stable isotope labeled compounds have an isotopic purity of 98 to 99%. Keeping the setting at 100% is appropriate for these compounds.
	Default: 100
2. Pattern Analysis	
Mass Tolerance [ppm]	Specifies the mass tolerance for the isotope search and XIC trace generation.
	Default: 5 ppm
Intensity Tolerance [%]	Specifies the relative intensity tolerance for isotope pattern matching.
	Default: 30
Intensity Threshold [%]	Specifies the isotope intensity threshold, relative to the pattern's base peak, for the theoretical isotope pattern.
	Default: 0.1
S/N Threshold	Specifies the signal-to-noise threshold for isotope pattern
S/IV Incshold	matching.

 Table 39.
 Analyze Labeled Compounds node (Sheet 2 of 3)

Parameter	Description
3. General Settings	
Mark Irregular Exchange	Adds the Irregular Exchange status to the following result table columns: Labeling Status column of the Compounds Table and the Status column of the Compounds per File Table.
	The Irregular Exchange status applies to compounds where the exchange rate for the mid-mass isotopologues is less than 5% of the exchange rate for the isotopologues.
	Default: True
Exclude Blanks	Specifies whether to exclude blank samples (Sample Type: Blank) from the main Compounds table.
	Default: True
Hide Unprocessed	Specifies whether to hide compounds with unassigned formulas in the Compounds table. The node does not process compounds without formulas.
	Default: True

 Table 39.
 Analyze Labeled Compounds node (Sheet 3 of 3)

Related Topics

• Working with the Isotopologues Distribution Chart

Detect Compounds Node

Use the Detect Compounds node to detect unknown compounds. The processing results for the Detect Compounds node appear in the Compounds per File table and the Features table.

For information about an untargeted processing workflow, see "Untargeted Workflows for Identifying Unknown Compounds" on page 24.

Table 40 describes the parameters for the Detect Compounds node.

 Table 40.
 Detect Compounds node parameters (Sheet 1 of 3)

Parameter	Description
1. General Settings	
Mass Tolerance	Specifies the mass tolerance for the XIC traces.
	Default: 5.0 ppm; range: 1–20.0 ppm
Intensity Tolerance	Specifies the intensity tolerance for the isotope pattern search.
	Default: 30%; range: 0 to 100%

Parameter	Description
S/N Threshold	Specifies the signal-to-noise threshold for mass spectral peaks (centroids). The unknown compound detection algorithm ignore centroids below this user-specified threshold.
	Default: 3; range: 0 to no upper limit
Min. Peak Intensity	Specifies the minimum base peak height to detect peaks in the XIC traces. The analysis does not report chromatographic peaks below this minimum peak intensity (peak height at the apex).
	The optimal minimum peak intensity setting depends upon the instrument (see Table 41).
	Default: 10 000
Ions	Specifies the adduct ions that might be in your samples.
	Default: [M+H] ⁺¹ , [M+K] ⁺¹ , [M+Na] ⁺¹
	Selection: The ion definitions in your Ion Definitions library
Base Ions	Specifies the adduct ions that you expect to have the highest intensity in your samples.
	Default: $[M+H]^{+1}$ and $[M-H]^{-1}$
	workflows (in the Common Workflows > Workflow Templates > E specify the following base ions: $[M+H]^{+1}$, $[M+NH_4]^{+1}$, and $[M-$
leachable compounds	ons includes the ammonium adduct because extractable and tend to be nonpolar compounds that require ammonium acetate as er for optimal chromatography.
leachable compounds a mobile phase modifi	tend to be nonpolar compounds that require ammonium acetate as
leachable compounds a mobile phase modifi	tend to be nonpolar compounds that require ammonium acetate as er for optimal chromatography. Specifies the minimum number of each possible element. The
leachable compounds a mobile phase modifi Min Element Counts	tend to be nonpolar compounds that require ammonium acetate as er for optimal chromatography. Specifies the minimum number of each possible element. The node uses these values for the isotope pattern search.
leachable compounds a mobile phase modifi Min Element Counts	tend to be nonpolar compounds that require ammonium acetate as er for optimal chromatography. Specifies the minimum number of each possible element. The node uses these values for the isotope pattern search. Default: C H Specifies the maximum number of each possible element. The
leachable compounds a mobile phase modifi Min Element Counts Max Element Counts	 tend to be nonpolar compounds that require ammonium acetate as er for optimal chromatography. Specifies the minimum number of each possible element. The node uses these values for the isotope pattern search. Default: C H Specifies the maximum number of each possible element. The node uses these values for the isotope pattern search.
leachable compounds	 tend to be nonpolar compounds that require ammonium acetate as er for optimal chromatography. Specifies the minimum number of each possible element. The node uses these values for the isotope pattern search. Default: C H Specifies the maximum number of each possible element. The node uses these values for the isotope pattern search.

Table 40. Detect Compounds node parameters (Sheet 2 of 3)

Parameter	Description
Max. Peak Width [min]	When Filter Peaks is set to True, specifies the maximum peak width at half height, in minutes, for detected chromatographic peaks.
	Default: 0.3; range: 0.05 to no upper limit
Remove Singlets	If set to True, does not report a component if the chromatographic peak for the component contains only A0 centroids.
	Default: True
Min. #Scans per Peak	Specifies the minimum number of scans (data points) required to define a chromatographic peak.
	Default: 5; range: 3 to 20
Min. #Isotopes	Specifies the minimum number of isotopes required in the mass spectral scans that define the chromatographic peak.
	Default: 1

Table 40. Detect Compounds node parameters (Sheet 3 of 3)

Table 41 lists the recommended range for the minimum peak intensity parameter. The optimal setting depends on the sensitivity of the mass spectrometer.

Table 41. Recommended minimum peak intensity range

Mass spectrometer	Minimum peak intensity (chromatographic peak height)
Q Exactive, Q Exactive Plus, Q Exactive HF	500 000 to 1 000 000
Orbitrap Fusion, Orbitrap Lumos, Orbitrap ID-X	50 000 to 100 000
Exactive, Exactive Plus, Orbitrap Elite, Orbitrap Velos Pro	100 000 to 500 000
LTQ Orbitrap XL, LTQ Orbitrap Velos	25 000 to 100 000

Fill Gaps Node

Use the Fill Gaps node to find chromatographic peaks that were detected by the Detect Compounds node in one of the input files but were missing from other input files in the file set.

For information about adding the Fill Gaps node to a processing workflow, see "Marking Background Compounds in an Untargeted Analysis" on page 28.

The Fill Gaps node adds the (hidden) Gap Status column to the Compounds Table and creates the related (hidden) Filled Gaps Table.

The Filled Gaps table describes how the node calculated the missing chromatographic peak areas. Clicking a row in the Filled Gaps table displays the gap-filled trace and the integrated The Fill Gaps node calculates the area of missing chromatographic peaks as follows.

- 1. Calculates the detection limit for each missing ion as an area of a simulated Gaussian peak that starts and ends at a zero intensity baseline. To calculate the area of the Gaussian peak, the node uses the expected peak width and the maximum spectrum noise in the expected retention time range multiplied by the S/N threshold.
- 2. Searches for the missing ion with the expected $m/z \times RT$ dimensions against all detected ions (in the mass list generated by the Detect Compounds node) while ignoring the assigned adduct type. If it finds a match (ion with the expected $m/z \times RT$ dimensions), it uses the ion's area to fill the gap and displays **Filled by Matching Ion** for the Fill Status.
- 3. If the node does not find a matching ion, it attempts to detect the peak at a lower intensity threshold using the Parameterless Peak Detection (PPD) algorithm. If it detects a chromatographic peak at a lower threshold, it uses the integrated peak area to fill the gap and displays **Re-detected Peak** for the Fill Status.
- 4. If the node does not find a chromatographic peak by using a lower intensity threshold, it fits a Gaussian peak to the XIC trace for the expected *m/z* range and displays **Filled by Simulated Peak** for the Fill Status.
- 5. If the filled area is still zero or lower than the detection limit, the node uses the detection limit value to fill the gap and displays **Filled by Spectrum Noise** for the Fill Status.
- 6. If the node cannot fill the gap, it displays Area Could Not Be Filled for the Fill Status.

Table 42 describes the parameters for the Fill Gaps node.

Table 42. Fill Gaps node parameters (Sheet 1 of 2)

Parameter	Description	
1. General Settings		
Mass Tolerance	Specifies the mass tolerance that the node uses to search for matching mass peaks.	
	Search range = expected mass ± mass tolerance/1e6	
	Default: 5 ppm; range: 0.1 to 20 ppm	
S/N Threshold	Specifies the minimum signal-to-noise threshold for centroids.	
	Default: 1.5; minimum: 1	

Parameter	Description
Use Real Peak Detection	Specifies whether the analysis fills gaps in the chromatogram by using a peak detection algorithm or only a peak simulation algorithm.
	True—The analysis uses the Parameterless Peak Detection (PPD) algorithm to fill the gaps with redetected low-intensity peaks. Using PPD can significantly increase the processing time.
	False—The analysis uses only a peak simulation algorithm to fill the gaps with simulated chromatographic peaks.

Table 42. Fill Gaps node parameters (Sheet 2 of 2)

Related Topics

• Filled Gaps Table

Group Compounds Node

Use the Group Compounds node to combine unknown compounds across the input file set by their molecular weight and retention time. This node also selects the best representative MS1scan and fragmentation tree, which the Predict Compositions node and search nodes use for identification.

Table 43 describes the parameters for the Group Compounds node.

Table 43. Group Compounds node parameters

Parameter	Description
1. Compound Consolidation	
Mass Tolerance	Specifies the mass tolerance that the node uses to search for matching mass peaks.
	Default: 5 ppm; range: 0.1 to 20 ppm
RT Tolerance [min]	Specifies the retention time tolerance, in minutes, that the node searches for mass peaks within the specified mass tolerance.
	Default: 0.05; range: 1.0
2. Fragment Data Select	tion
Preferred Ions	Select the preferred ions (adducts) from the list. The application uses the list to select the best fragmentation data for each compound to submit to an mzCloud or mzVault search.
	Selections: One or more of the ions in your Ion Definitions library
	Default: [M+H]+1 and [M–H]–1

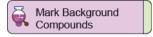
Peak Area Refinement Nodes

Use these nodes to hide background compounds or to normalize the chromatographic peak areas:

- Mark Background Compounds Node–Unknown Compounds
- Normalize Areas Node

Mark Background Compounds Node–Unknown Compounds

Use the Mark Background Compounds node to identify the background compounds, which are compounds also found in the blank samples, in the Compounds Table.



For information about the parameter settings, see "Mark Background Compounds Node-Expected Compounds" on page 240.

Always place the Mark Background Compounds node at the end of the processing workflow by following these examples:

- Group Compounds > Mark Background Compounds
- Group Compounds > Fill Gaps > Mark Background Compounds
- Group Compounds > Fill Gaps > Normalize Areas > Mark Background Compounds

IMPORTANT As you drag the Fill Gaps and the Normalize Areas nodes into the Workflow Tree pane, the application automatically connects the Group Compounds node to the Fill Gaps node, and the Fill Gaps node to the Normalize Areas node. The application does not connect the Mark Background Compounds node to other nodes, so you must manually make the appropriate connections. For more information about the node connections, see "Marking Background Compounds in an Untargeted Analysis" on page 28.

Normalize Areas Node

Use the Normalize Areas node to normalize the chromatographic peak areas when you need the analysis to compensate for batch effects.

When the Normalize Areas node is part of a processing workflow, the Use Normalized Areas check box is available in the Descriptive Statistics view and the Principal Component Analysis view. The Normalize Areas node adds the Normalized Area column to the Compounds Table. When the analysis includes quality control samples, the node also adds the "QC" columns in the Compounds table. You can view the corrected compound areas in the Compound Area Corrections view (see "Viewing the QC-Based Compound Area Correction Curves" on page 367).

IMPORTANT When the Normalize Areas and Mark Background Compounds nodes are part of a processing workflow, connect the Normalize Areas node to the Mark Background Compounds node.

If the input files include Quality Control samples, the application automatically performs a batch normalization on each compound that meets the user-specified criteria.

Table 44 describes the parameters for the Normalize Areas node.

Table 44. Normalize Areas node parameters (Sheet 1 of 3)

Parameter	Description	
	-	

1. QC-based Area Correction

The analysis uses the QC samples to create a regression curve for each detected compound. The analysis does not create a regression curve for a particular compound or does not correct the areas in the non-QC samples unless all three of these conditions are met:

- It detects the compound in the specified minimum percentage of QC samples.
- The relative standard deviation of the detected peak areas for the compound in the QC samples does not exceed the specified threshold.
- The number of samples acquired between the QC samples does not exceed the specified number.

If the analysis does not create a regression curve for a compound, it does not perform a batch normalization of the peak areas for the compound in the non-QC samples (Sample type: Sample, Control, or Standard).

Regression Model	Specifies the regression model as Linear or Cubic Spline.
	Default: Linear
Min. QC Coverage [%]	Specifies the minimum percentage of the QC samples where the analysis must detect a particular compound before it creates a regression curve for the compound. If the coverage falls below this value, the analysis does NOT perform a batch normalization of the peak areas for the compound in the non-QC samples. Default: 50%; range: 25 to 100%

Parameter	Description
Max. QC Area RSD [%]	Specifies the maximum relative standard deviation (RSD%) for the areas for a particular compound across the QC samples. If the RSD% exceeds this maximum percentage, the analysis does not create a regression curve for the compound.
	The application checks the RSD% of the peak areas both before and after applying a correction for the relative acquisition time. If the RSD% for the corrected peak areas differs by more than 75% of the user-specified threshold, the application does not create a corrected-area regression curve for the compound.
	Default: 30%; range: 10 to 50%
Max. #Files Between QC Files	The application checks the acquisition time stamp for each input file in an analysis and orders the samples within an analysi set by acquisition time.
	Specifies the maximum number of non-QC samples that you can acquire between the QC samples. If the application detects more non-QC samples between the QC samples than the maximum allowable number, it does not correct the chromatographic peak areas in the non-QC samples—that is, it does not perform a batch normalization on these samples.
	Default: 15; range: 1 to unchecked value
2. Area Normalization	
Normalization Type	For selections other than None, the analysis normalizes the chromatographic peaks areas after performing a batch normalization based on the QC samples.
	 Selections: None (Default)—The analysis does not normalize the chromatographic peak areas for injection variability. Constant Sum Constant Mean Constant Median Median Absolute Deviation (MAD) Scaling Factor—The node multiplies the peak areas by the numeric values for the user-specified study factor.
	Tip For best results, select Constant Mean for input file sets that include matrix blanks.

Table 44.	Normalize Areas node parameters (Sheet 2 of 3)
-----------	--

Parameter	Description
Exclude Blanks	Specifies whether to exclude Blank sample types from normalization.
	Default: True
	Tip For best results, follow these guidelines:
	• Select True for solvent blanks to exclude them from the normalization process. Otherwise, the Fill Gaps node adds small "noise" peaks to the solvent blanks for each detected compound in the input file set. The Normalize Areas node magnifies these small peaks, causing the sample-to-blank ratio to fall below the user-specified value in the Mark Background Compounds node. The Mark Background Compounds node then hides most of the detected compounds across the input file set (marks them as background compounds).
	• Select False for matrix blanks, such as plasma and urine, as these blanks typically contain a large number of compounds that you might want to hide.
3. Scaling Factor	
Study Factor Name	Specifies the study factor to use as the scaling factor for the chromatographic peak areas when you select Scaling Factor in the Normalization Type list.
	Multiplies a compound's peak area by the input file's value for the selected numerical study factor.
	Selections: Defined numerical study factors Default: Empty

Table 44.	Normalize Areas node parameters (Sheet 3 of 3)
-----------	--

Compound Identification Nodes

Use these nodes to identify unknown compounds:

- Assign Compound Annotations Node
- Predict Compositions Node
- Search ChemSpider Node
- Search Mass Lists Node
- Search mzCloud Node
- Search mzVault Node

Assign Compound Annotations Node

Use the Assign Compound Annotations node to select the preferred data sources for the following annotations: name, formula, and structure. The application attempts to assign the annotations provided by the first data source. If the first source does not provide the annotation, the application uses the second data source, and so on. If the processing workflow does not include the Assign Compound Annotations node, the application does not populate the Name, Formula, or Structure columns of the Compounds table.

Table 45 describes the parameters for the Assign Compound Annotations node.

Table 45.	Assign Compound	Annotation node	parameters (Sheet 1 of 2)

Parameter	Description
1. General Settings	
Mass Tolerance	Specifies the mass tolerance for validating the annotations.
	Default: 5 ppm; range: 0.1–20 ppm

2. Data Sources

Available sources: mzCloud Search, mzVault Search, BioCyc Search, Mass List Match, ChemSpider Search, and Predicted Compositions

Data Source #1	Specifies the primary source for the compound annotations.	
	Default: mzCloud Search	
	Tip If a stable isotope analysis misidentifies known compounds, consider reprocessing the analysis after selecting your custom mass list as the first data source.	
Data Source #2	Specifies the secondary source for compound annotations if the primary source is unavailable.	
	Default: Predicted Compositions	

Parameter	Description
Data Source #3	Specifies the source for compound annotations if the primary and secondary sources are unavailable.
	Default: Mass List Search
Data Source #4	Specifies the source for compound annotations if all the previous sources are unavailable.
	Default: ChemSpider Search
Data Source #5–7	Use to specify additional annotation sources in the list.
	Default: Empty
3. Scoring Rules	
Use mzLogic	When set to True, uses the score from the Apply mzLogic node to select the best candidate.
	Default: True
Use Spectral Distance	When set to True, uses the SFit score from the Apply Spectral Distance node to select the best candidate.
	Uses the SFit Threshold and SFit Range values to limit the number of valid candidates.
	Default: True
SFit Threshold	Specifies the minimum SFit score for a candidate.
	Default: 20; range: 0 to 100
SFit Range	Specifies the maximum allowed difference between the SFit scores for the best and worst candidates.
	Default: 20; range: 0 to 100

 Table 45.
 Assign Compound Annotation node parameters (Sheet 2 of 2)

Predict Compositions Node

Use the Predict Compositions node to predict the chemical formulas of the unknown compounds. This node creates the Predicted Compositions Table and populates the Formula column of the Compounds Table when the processing workflow also includes the Assign Compound Annotations node.

Table 46 describes the parameters for the Predict Compositions node.

Parameter	Description	
1. Prediction Settings		
Mass Tolerance	Specifies the mass tolerance for the XIC traces.	
	Default: 5.0 ppm; range: 1–20.0 ppm	
Min. Element Counts	Specifies the minimum count for each element in the hypothetical compound. If an element is not listed, its minimum count is zero.	
	Default: C H	
Max. Element Counts	Specifies the maximum count for each element in the hypothetical compound. If an element is not listed, its maximum count is the same as its minimum count.	
	Default: C90 H190 Br3 Cl4 K2 N10 Na2 O15 P2 S5	

Table 46. Predict Compositions node parameters (Sheet 1 of 4)

Table 46.	Predict Compositions	node parameters	(Sheet 2 of 4)
	i iouior oonipooniono	nouo paramotoro	

Parameter	Description
Min. RDBE and Max. RDBE	Specifies a range of values for ring and double-bond equivalents. The RDBE value is a measure of the number of unsaturated bonds in a compound. The specified value limits the calculated formulas to only those that make sense chemically.

The following formula determines the RDBE value for an elemental composition:

$$D = 1 + \frac{\left[\sum_{i=1}^{i\max} Ni(Vi-2)\right]}{2}$$

where:

	• <i>D</i> is the value for the RDB equivalents
	• <i>i</i> max is the total number of different elements in the composition
	• <i>Ni</i> is the number of atoms of element i
	• <i>Vi</i> is the valence of atom i
Min. H/C	Specifies the minimum hydrogen-to-carbon ratio.
	Default: 0.1
	The value of 0 means no limit. The application does not accept negative values.
Max. H/C	Specifies the maximum hydrogen-to-carbon ratio.
	Default: 3
	The value of 0 means no limit. The application does not accept negative values.

Parameter	Description
Note For most compout to 2.0.	nds, the hydrogen-to-carbon ratio falls within the range from 0.5
• Long chain alkanes	have an H/C ratio of approximately 2.
Polycyclic aromatics	have an H/C ratio of approximately 0.5.
Max. #Candidates	Specifies the maximum number of compositions to store for each compound in the result file.
	Default: 10; range: 1–50
Max # Internal Candidates	Specifies the maximum number of compositions to calculate for each detected component.
	Default: 200; range: 1 to unchecked
	Tip To optimize the processing time, the maximum number of internal candidates is limited to 200 by default and sorted by the mass error. Normally, this works well for masses below 500 Da. However, as a component's mass or the number of elements increases, the number of possible elemental compositions also increases. This increase means that the chance of rejecting the correct formula, when it has a higher mass error than the first 200 candidates, also increases.
	For samples with components that have a relatively large number of elements or a mass above 500 Da, consider increasing the limit to 500.
2. Pattern Matching	
Intensity Tolerance [%]	Specifies the intensity tolerance for the isotope pattern search.
	Default: 30%; range: 0 to 100%
Intensity Threshold [%]	Specifies the intensity threshold, relative to the base peak (most intense ion) in the isotope pattern, for the isotope pattern search The analysis ignores isotopes below this threshold.
	Default: 0.1%; range: 0.1 to 10%
S/N Threshold	Specifies the signal-to-noise threshold for the isotope search. Isotopes with a theoretical intensity below the threshold are not required.
Min. Spectral Fit [%]	Specifies the minimum spectral fit for reporting a predicted composition in the Predicted Compositions result table.
	composition in the redicted compositions result table.

Table 46. Predict Compositions node parameters (Sheet 3 of 4)

Parameter	Description
Min. Pattern Cov. [%]	Specifies the minimum percentage for the summed intensity of the matching isotope peaks in the measured MS1 spectrum relative to the summed intensity of the theoretical isotope pattern.
	$\frac{Summed \ intensity \ of \ the \ matching \ isotope \ peaks \times 100}{Summed \ intensity \ of \ the \ theoretical \ isotope \ pattern}$
Use Dynamic Recalibration	Specifies whether the application uses the dynamic recalibration algorithm to shift the theoretical pattern for the candidate formula by the difference in the observed m/z value of the leftmost (A0) isotopic peak in the measured spectrum.
	Use dynamic recalibration when there is a systematic error (due to calibration) in the measured spectrum.
	Default: True
3. Fragments Matching	
Use Fragments Matching	Specifies whether the application uses the fragment matching algorithm, which ranks the identified candidates (chemical formulas) by the number of matching centroids (with an m/z value that matches a subset of the elemental composition for a particular candidate) in the fragmentation scan for the precursor ion. Default: True
Mass Tolerance	Specifies the mass tolerance for matching the centroids in the
	fragmentation scans to the <i>m/z</i> values for the expected fragments Default: 5 ppm; range: 0 to unchecked
S/N Threshold	Specifies the signal-to-noise threshold for the centroids in the fragmentation scans. The node ignores centroids with an intensity below this threshold.
	Default: 3; range: 0 to unchecked

 Table 46.
 Predict Compositions node parameters (Sheet 4 of 4)

Search ChemSpider Node

Use the Search ChemSpider node to search mass spectral databases for matching compounds within a specified mass tolerance range or with a certain elemental composition. This node requires input from the Group Compounds node. It adds the #ChemSpider Results column to the Compounds Table, and creates the ChemSpider Results Table.

When the processing workflow includes a ChemSpider search, the processing computer must have Internet access. To verify whether the processing computer can access the ChemSpider database, run the Communication test. See Chapter 17, "Testing Communication to the Online Databases."

Table 47 describes the parameters for the Search ChemSpider node.

Parameter	Description
1. Search Settings	
Database(s)	Specifies the databases for the ChemSpider search.
	Default: KEGG
Search Mode	Default: By Formula or Mass
	Selections: By Formula or Mass, By Formula and Mass, By Formula or Mass, By Mass Only
Mass Tolerance	Specifies the mass tolerance that the node uses to search for matching mass peaks.
	Default: 5 ppm; range: 0.1 to 20 ppm or 0.0 to 0.1 Da
Max# of Results Per Compound	Specifies the maximum number of hits (matches) to return (store in the result file).
	Default: 100; range: 1 to 2000
Max # of Predicted Compositions to be	Specifies the maximum number of predicted compositions to search for in the ChemSpider database.
Searched per Compound	Default: 3; range: 1 to 100

Table 47. Search ChemSpider node parameters (Sheet 1 of 2)

Parameter	Description	
Result Order (for Max#	Specifies the sort order in the ChemSpider Results table.	
of Results per Compound)	Selections:	
	• Order By Reference Count (DESC)—Sorts the search results by the number of references for each compound.	
	• Order By Data Source Count (DESC)—Sorts the search results by the number of data sources.	
	• Order By Mass Deviation (ASQ)—Sorts the search results by mass deviation from the expected mass.	
	• Order By PubMed Count (DESC)—Sorts the search results by the number of PubMed references.	
	• Order By RSC Count (DESC)—Sorts the search results by the number of RSC references.	
	• Order By CSID (ASQ)—Sorts the search results by the ChemSpider ID.	
2. Predicted Composition	Annotation	
Check All Predicted Compositions	Specifies whether to add a flag to the Predicted Compositions table. When set to True, the Search ChemSpider node adds the Ir ChemSpider column to the Predicted Compositions table and marks the matched Predicted Compositions with an X.	
	Default: False	

Table 47.	Search ChemS	pider node	parameters (Sheet 2 of 2)
-----------	--------------	------------	---------------------------

Search Mass Lists Node

Use the Search Mass Lists node to search mass lists for masses that match the detected compounds. This node requires input from the Group Compounds node, adds the #Matched Masses to the Compounds Table, and creates the Mass List Search Results Table.

Table 48 describes the parameters for the Mass List Search node.

Parameter	Description	
1. Search Settings		
Mass Lists	Specifies the mass list files that the node searches for matching masses.	
	The mass list must have the following columns—Mass, Retentio Time, and Name. The mass list can also contain the following additional columns—Molecular Structure and Text Annotation.	
	 To select input files for the mass list search 	
	1. In the Mass box, click the browse icon,	
	The Select Input Files dialog box opens.	
	2. Select the check boxes for the files that you want to use for the mass list search.	
	3. Click OK .	
Use Retention Time	Specifies whether to search for compounds by retention time in addition to mass.	
	Default: True	
RT Tolerance [min]	Specifies the retention time tolerance, in minutes, for the search. The node searches for matching peaks in a retention time window equal to the expected retention time plus or minus the specified RT tolerance.	
	Default: 0.05; range: 1.0	
Mass Tolerance	Specifies the mass tolerance that the node uses to search for matching compounds.	
	Default: 5 ppm; range: 0.1 to 20 ppm or 0.0 to 0.1 Da	

 Table 48.
 Search Mass List node parameters

Table 49 describes the table columns in the Select Input Files dialog box.

Table column	Description
Selected	Selecting the check box for a mass list adds the list to the search.
Filename	Displays the file name of the imported file.
Description	Editable text field.
	Clicking the table cell displays a text entry field for typing a name or a description of the mass list. Use this column to name and sort your mass lists.
File Size	Displays the file size of the imported file.
Uploaded	Displays the date (month/day/year) and time (hour/minute) when you added the file to the library in the following format:
	MM/DD/yyyy HH:mm
Updated	Displays the date and time when the file was updated.
Context	Displays the source of the mass list—for example, Import from CSV or Import from XML.
State	Specifies whether the mass list is available, corrupted, or missing.
	If you remove a mass list from the ServerFiles folder or edit a mass list in the ServerFiles Folder, and then restart the application, the mass list's state changes to Missing or Corrupted, respectively.

Table 49. Select Input File(s) dialog box parameters

Search mzCloud Node

Use the Search mzCloud node to search the mzCloud database for matching fragmentation spectra. This node creates the mzCloud Results Table, and adds the #mzCloud Results and mzCloud Best Match columns to the Compounds table.

The Search mzCloud node requires input from the Group Compounds node.

Note In addition to running an automated search with an analysis, you can manually submit a fragmentation scan to the mzCloud database from the Mass Spectrum view for an active result file.

When the processing workflow includes an mzCloud search, the processing computer must have Internet access. To verify whether the processing computer can access the mzCloud database, run the Communication tests.

Table 50 describes the parameters for the Search mzCloud node.

Parameter	Description
1. General Settings	
Compound Classes	Specifies the compound classes for the search. Select All or select specific classes.
	Default: All
Precursor Mass Tolerance	Specifies the tolerance for the precursor mass.
	Default: 10 ppm; range: 0–0.1 Da or 0–20 ppm
FT Fragment Mass Tolerance	Specifies the mass tolerance for high-resolution fragmentation scans performed in the Orbitrap analyzer (FTMS).
	Default: 10 ppm; range: 0–0.5 Da or 0–100 ppm
IT Fragment Mass Tolerance	Specifies the mass tolerance for low-resolution fragmentation scans performed in the ion trap analyzer (ITMS).
	Default: 0.4 Da; range: 0–0.1 Da or 0–20 ppm
Library	Specifies the mzCloud library for the search: Autoprocessed, Reference, or both.
	• Reference library—Contains only spectra that have been manually curated by mass spectrometry experts
	• Autoprocessed library—Contains spectra curated with an automated process. As time permits, spectra in the Autoprocessed library are manually curated and transferred to the Reference library.
Post Processing	Specifies whether to search the library for filtered or recalibrated spectra.
	Selections:
	• Filtered—Removes extraneous mass peaks that do not match the theoretical mass spectrum.
	• Recalibrated (Default)—The mass peaks for the known ion fragments are recalibrated to match the theoretical mass spectrum by a series of manually supervised ion calibration steps.

Table 50. Search mzCloud node parameters (Sheet 1 of 5)

Parameter	Description
Max. # Results	Specifies the maximum number of hits for each compound and search spectrum to store in the result file
	Default: 10; range: 1 to 50
Annotate Matching Fragments	Specifies whether the processing workflow annotates the matching fragments with structures that you can review in the Spectrum view of the result page.
	Default: False
2. DDA Search	
Identity Search	Specifies the identity search algorithm.
	Selections:
	• Confidence—Takes into account the dot product of both spectra, distribution of peaks, activation energy difference, and polarity, and uses a machine-learning Bayesian Network model to estimate the likelihood of a correct match.
	 HighChem DP—Alternative score using a modified dot product algorithm that includes a "hard penalty" function for spectra with a low number of fragment ions.
	• HighChem HighRes—Use for high-resolution data. Correlation algorithm that uses the geometric mean of a modified Spearman's rank order correlation to separately determine the intensity and <i>m/z</i> accuracy of the fragment ions.
	• NIST
	• Cosine—Dot product algorithm with fragment intensities weighted by 0.75 and no weighting on fragment <i>m/z</i> values.
	Default: HighChem HighRes
Match Activation Type	Specifies whether to only search for fragmentation spectra with the same activation type.
	Default: True

Table 50. Search mzCloud node parameters (Sheet 2 of 5)

Parameter	Description
Match Activation Energy	Specifies whether to search for fragmentation spectra generated by the same ion activation energy within a tolerance or by any ion activation energy.
	Default: Match with Tolerance
Activation Energy Tolerance	Specifies the tolerance as an absolute value for the ion activation energy used to generate the fragmentation spectrum.
	For example, if the ion activation energy used to generate your spectrum was a normalized collision energy of 35% and you specify an ion activation energy tolerance of ± 20 the search looks for spectra with an ion activation energy from 15 to 55.
	Default: 20; range: 0 to 200
Apply Intensity Threshold	Specifies whether to apply an automatic intensity threshold that sets the threshold intensity by calculating the spectrum noise level.
	Default: True

Table 50. Search mzCloud node parameters (Sheet 3 of 5)

Parameter	Description
Similarity Search	Specifies the similarity search algorithm.
	Selections:
	• Confidence Forward— Takes into account the dot product of the forward search match and distribution of peaks, and the ratio of the most intensive matching peaks. A higher similarity score indicates the extent to which the unknown component resembles the library compound.
	• Confidence Reverse—Takes into account the dot product of the reverse search match and distribution of peaks, and the ratio of the most intensive matching peaks. A higher similarity score indicates the extent to which the library compound resembles the unknown component.
	• None (Default)—Does not run a similarity search.
	• Similarity Forward—Searches for a match between the best fragmentation scan for a compound (across the input file set) and a fragmentation scan in the mzCloud database. Unlike the Identity Search, this search ignores the <i>m</i> / <i>z</i> value of the precursor ion.
	• Similarity Reverse—Searches for a match between th fragmentation scans in the mzCloud database and th best fragmentation scan for a compound (across the input file set).
Match Factor Threshold	Specifies the minimum match factor for reporting a spectrum match.
	Default: 60; range: 0 to 100%
3. DIA Search	
Use DIA Scans for Search	Specifies whether to submit data independent scans to th mzCloud database for a spectral search.
	Default: False
Max. Isolation Width [Da]	Specifies the maximum MS2 isolation width. The isolation width for a scan is listed in its scan header.
	Default: 500 Da; minimum width: 10 Da

Table 50.	Search mzCloud	node parameters	(Sheet 4 of 5)
-----------	----------------	-----------------	----------------

Parameter	Description
Match Activation Type	See Match Activation Type under DDA Search.
	Default: False
Match Activation Energy	See Match Activation Energy under DDA Search.
	Default: Any
Activation Energy Tolerance	See Activation Energy Tolerance under DDA Search.
	Default: 100
Apply Intensity Threshold	See Apply Intensity Threshold under DDA Search.
	Default: False
Match Factor Threshold	See Match Factor Threshold under DDA Search.
	Default: 20

Table 50. Search mzCloud node parameters (Sheet 5 of 5)

Related Topics

• mzCloud Results Table

Search mzVault Node

Use the Search mzVault node to search a local mass spectra database for compounds of interest. This node requires input from the Group Compounds node and creates the mzVault Results table.

To use mzVault libraries created with the mzVault 1.1 or earlier application

Change the following settings:

- For the Match Ion Activation Type, select **False**.
- For the Match Ion Activation Energy, select **Any**.
- For the Match Ionization Method, select False.

Table 51 describes the parameters for the Search mzVault node.

Table 51. Search mzVault node parameters (Sheet 1 of 4)

Parameter	Description
1. Search Settings	
mzVault Library	Specifies the mzVault libraries to search. You can select one or more of the mzVault libraries in your Spectral Libraries list.

Parameter	Description
Compound Classes	Specifies the compound classes for the search. Select All or select specific classes.
	Default: All
Match Ion Activation Type	Specifies whether to only search for library scans that match the ion activation type of the query spectrum.
	Default: True
Match Ion Activation Energy	Specifies whether to search for library scans generated by the same ion activation energy within a tolerance of the query spectrum or by any ion activation energy.
	Default: Match with Tolerance
Ion Activation Energy Tolerance	Specifies the tolerance as an absolute value for the ion activation energy used to generate the fragmentation spectrum.
	For example, if the ion activation energy used to generate your spectrum was a normalized collision energy of 35% and you specify an ion activation energy tolerance of ± 20 , the search looks for spectra with an ion activation energy from 15 to 55.
	Default: 20; range: 0 to 200
Match Ionization Method	Specifies whether to only search for library spectra from the same ionization method (for example, ESI, APCI, and so on) as the query spectrum.
	Default: True
Apply Intensity Threshold	Specifies whether to apply an automatic intensity threshold that sets the threshold intensity by calculating the spectrum noise level.
	Default: True
Remove Precursor Ion	When set to True, the search ignores (removes from consideration) mass peaks within 2.2 Da of the precursor ion's m/z value in the query spectrum—that is, the match score is not negatively affected if the library spectrum
	does not include a mass peak within 2.2 Da of the m/z value for the precursor ion.

Table 51. Search mzVault node parameters (Sheet 2 of 4)

Parameter	Description
Precursor Mass Tolerance	Specifies the tolerance for the precursor mass.
	Default: 10 ppm; range: 0–0.1 Da or 0–20 ppm
FT Fragment Mass Tolerance	Specifies the mass tolerance for high-resolution fragmentation scans performed in the Orbitrap analyzer (FTMS).
	Default: 10 ppm; range: 0–0.5 Da or 0–100 ppm
IT Fragment Mass Tolerance	Specifies the mass tolerance for low-resolution fragmentation scans performed in the ion trap analyzer (ITMS).
	Default: 0.4 Da; range: 0–0.1 Da or 0–20 ppm
Match Analyzer Type	Specifies whether to only search for library spectra from the same mass analyzer type as the query spectrum.
	Default: True
Search Algorithm	Specifies the identity search algorithm.
	Selections:
	 HighChem DP—Alternative score using a modified dot product algorithm that includes a "hard penalty function for spectra with a low number of fragment ions.
	• HighChem HighRes—Use for high-resolution data Correlation algorithm that uses the geometric mean of a modified Spearman's rank order correlation to separately determine the intensity and <i>m/z</i> accuracy of the fragment ions.
	• NIST
	Default: HighChem HighRes
Match Factor Threshold	Specifies the minimum match factor for reporting a spectrum match.
	Default: 50; range: 0 to 100%

Table 51. Search mzVault node parameters (Sheet 3 of 4)

Parameter	Description
Max. # Results	Specifies the maximum number of hits for each compound to store in the result file. Displays the hits with the highest match score above the cutoff storage number in the result file.
	Default: 10
RT Tolerance [min]	Specifies the retention time tolerance for the library search.
	Default: 2; range: 0 to 10 minutes
Use Retention Time	Specifies whether to filter the database hits by their retention time. Filters out scans without retention time information.
	Default: False

Table 51. Search mzVault node parameters (Sheet 4 of 4)

Related Topics

• Modifying the Spectral Libraries List

Pathway Mapping Nodes

Use these nodes to map detected compounds to a biochemical pathway:

- Map to BioCyc Pathways Node
- Map to KEGG Pathways Node
- Map to Metabolika Pathways Node

Map to BioCyc Pathways Node

Use the Map to BioCyc Pathways node to map the BioCyc pathways for each compound. This node requires input from the Group Compounds node—that is, it requires a list of molecular weights or chemical formulas.

The Map to BioCyc Pathways node adds the following items to the result file:

- The BioCyc Pathways and BioCyc Results main tables
- The #BioCyc Pathways and BioCyc Pathways columns in the main Compounds table
- The BioCyc Compound IDs, BioCyc Compound Names, and BioCyc Compound Formula columns in the related Compounds table and the related BioCyc Pathways table.

Table 53 describes the parameters for the Map to BioCyc Pathways node.

Table 52.	Map to BioCyc	Pathways node	parameters
-----------	---------------	---------------	------------

Parameter	Description
1. Search Settings	
BioCyc Database/Organism to be Searched	Specifies the databases for the search.
Search Mode	Default: By Formula or Mass
	Selections: By Formula or Mass, By Formula and Mass, By Formula or Mass, By Mass Only
2. By Mass Search Settings	
Mass Tolerance	Specifies the mass tolerance that the node uses to search for matching compounds.
	Default: 5 ppm; range: 0.1 to 20 ppm
3. By Formula Search Settings	
Max. # of Predicted Compositions to be	Specifies the maximum number of predicted compositions to search for per compound.
Searched per Compound	Default: 3; range: 1.0
4. Display Settings	
Maximum # of Pathways in "Pathways" Column	Specifies the maximum number (n) of pathways to display by name in the Pathways column of the Compounds result table. The Pathways column displays the names of the first $n - 1$ pathways that include the compound. The other pathways that include the compound are grouped in the Other category.
	Default: 20; range: 1–30

Map to KEGG Pathways Node

The KEGG (Kyoto Encyclopedia of Genes and Genomes) pathway database contains connection diagrams of molecular interactions, reactions, and relations. Use the Map to KEGG Pathways node to add explanations to the result file about the reaction pathways for each detected compound. The node requires input from the Group Compounds node—that is, it requires a list of molecular weights or chemical formulas.

The Map to KEGG Pathways node adds the following items to the result file:

- The KEGG Pathways table
- The KEGG Compound IDs, KEGG Compound Names, and KEGG Compound Formula columns in the related Compounds Table
- The #Pathways and Pathways columns in the main Compounds table

Table 53 describes the parameters for the Map to KEGG Pathways node.

When the processing workflow includes the Map to KEGG Pathways node, the processing computer must have Internet access. To verify whether the processing computer can access the KEGG Pathways database, run the Communication test. See Chapter 17, "Testing Communication to the Online Databases."

Parameter	Description
1. Search Settings	
Search Mode	Specifies the search mode.
	Default: By Formula or Mass
	Selections: By Formula or Mass, By Formula and Mass, By Formula or Mass, By Mass Only
2. By Mass Search Settings	
Mass Tolerance	Specifies the mass tolerance that the node uses to search for matching compounds.
	Default: 5 ppm; range: 0.1 to 20 ppm
3. By Formula Search Settings	
Max. # of Predicted Compositions to be	Specifies the maximum number of predicted compositions to search for per compound.
Searched per Compound	Default: 3; range: 1.0

Table 53. Map to KEGG Pathways node parameters (Sheet 1 of 2)

Parameter	Description
4. Display Settings	
Maximum # of Pathways in "Pathways" Column	Specifies the maximum number (n) of pathways to display by name in the Pathways column of the Compounds result table. The Pathways column displays the names of the first $n - 1$ pathways that include the compound. The other pathways that include the compound are grouped in the Other category.
	Default: 20; range: 1–30

Table 53. Map to KEGG Pathways node parameters (Sheet 2 of 2)

Map to Metabolika Pathways Node

To search your local database of Metabolika pathways for pathways that include matching structures for the unknown compounds in your data set (by formula, mass, or both), add the Metabolika Pathways node to the processing workflow and select the pathways to search.

Note 378 Metabolika pathway files are automatically installed with the application. You can edit these pathways or create your own pathways by using the Metabolika pathway editor ("Modifying the Metabolika Pathways List" on page 528).

Table 54 describes the parameters for the Map to Metabolika Pathways node.

Parameter	Description
1. Search Settings	
Metabolika Pathways	Specifies the Metabolika pathways to search.
	Default: All
Search Mode	Specifies whether to search by formula, mass, or both.
	Default: By Formula or Mass
	Selections: By Formula and Mass, By Formula Only, By Formula or Mass, By Mass Only
2. By Mass Search Settings	
Mass Tolerance	Specifies the mass tolerance that the node uses to search for matching compounds.
	Default: 5 ppm; range: 0.1 to 20 ppm

Table 54. Map to Metabolika Pathways node parameters (Sheet 1 of 2)

Parameter	Description
3. By Formula Search Settings	
Max. # of Predicted Compositions to be Searched per Compound	Specifies the maximum number of predicted compositions to search for per compound. Default: 3; range: 1 to 100
4. Display Settings	
Maximum # of Pathways in "Pathways" Column	Specifies the maximum number (n) of pathways to display by name in the Pathways column of the Compounds result table. The Pathways column displays the names of the first n-1 pathways that include the compound. The other pathways that include the compound are grouped in the Other category. Default: 20; range: 1 to 30

Table 54. Map to Metabolika Pathways node parameters (Sheet 2 of 2)

Compound Scoring Nodes

Use these nodes to score the explanations for each detected compound.

- Apply Spectral Distance Node
- Apply mzLogic Node
- Calculate Mass Defect Node
- Compound Class Scoring Node
- Generate Molecular Networks Node (Beta)
- Pattern Scoring Node

To specify the isotope pattern for the Pattern Scoring node or the Pattern Trace node, see this topic "Setting Up Isotope Patterns for Pattern Scoring or Pattern Traces" on page 142.

Apply Spectral Distance Node

Use the Apply Spectral Distance node to calculate a spectral similarity score (SFit[%]) for the compound annotations. The SFit[%] column appears in the related "compound annotations" tables for the compound selected in the Compounds table. Depending on the processing workflow, the related "compound annotations" tables can include any of these tables: ChemSpider Results, Mass List Search Results, Metabolika Results, and BioCyc Results. In addition to providing an SFit score for you to review in the related "compound annotations" tables, the node sends the scores to the Assign Compound Annotation node.

Table 55 describes the parameters for the Apply Spectral Distance node.

Parameter	Description
1. General Settings	
Mass Tolerance	Specifies the mass tolerance that the node uses to search for matching mass peaks.
	Search range = expected mass ± mass tolerance/1e6
	Default: 5 ppm; range: 0.1 to 20 ppm
Intensity Tolerance [%]	Specifies the relative intensity tolerance of the mass spectral peaks in the isotope pattern.
	Default: 30; range: 0 to 100%
Intensity Threshold [%]	Specifies the isotope intensity threshold, relative to the base peak of the isotope pattern, that the node uses for pattern simulation. The node does not add isotopes below this threshold to the simulated pattern.
	Default: 0.1
S/N Threshold	Specifies the signal-to-noise threshold for the isotope search. The node does not include isotopes that are expected to be below this threshold in the SFit score—that is, these isotopes are not required in the measured isotope pattern.
	Default: 3
Use Dynamic Recalibration	Specifies whether to shift the theoretical isotope pattern if the pattern base peak in the query spectrum is shifted.
	Default: True

Table 55. Apply Spectral Distance node parameters

Related Topics

• ChemSpider Results Table

- Mass List Search Results Table
- BioCyc Results Table
- Metabolika Results Table

Apply mzLogic Node

Use the Apply mzLogic node to score explanations from the ChemSpider node, Search Mass List node, Map to BioCyc Pathways node, and Map to Metabolika Pathways node.

Table 56. Apply mzLogic node parameters

Parameter	Description
1. Search Settings	
Max # Compounds	Specifies the maximum number of compounds to display and score in the result table.
	Default: 0 (no maximum limit)
Max # mzCloud Similar Results to Consider per Compound	Specifies the maximum number of compounds to consider from an mzCloud similarity search. Increasing the number of compounds to consider increases the processing time.
	Range: 5 to 100
Match Factor Threshold	Specifies the minimum match score returned for a compound by an mzCloud similarity search. The analysis ignores compounds with match scores below this threshold.
	Default: 30
Advanced parameters	
FT Fragment Mass Tolerance	Specifies the mass tolerance for the mass peaks in high-resolution fragmentation spectra when searching the mzCloud spectral database.
	Default: 10 ppm; range: 0 to 0.5 Da or 0 to 100 ppm
IT Fragment Mass Tolerance	Specifies the mass tolerance for the mass peaks in low-resolution fragmentation spectra when searching the mzCloud spectral database.
	Default: 0.4 Da; range: 0 to 1 Da or 0 to 1500 ppm

Calculate Mass Defect Node

Use this node to calculate the mass defect for each detected compound. You can specify up to four calculation methods and up to five Kendrick formulas. This node adds the Mass Defect column to the Compounds table for reviewing the result of the specified calculations.

Table 31 describes the parameters for the Filter By Mass Defect node.

 Table 57.
 Filter By Mass Defect node parameters

Parameter	Description
1. Mass Defect	
Fractional Mass	Fractional Mass = exact mass - floor (exact mass)
Standard Mass Defect	Standard Mass Defect = exact mass – nominal mass
Relative Mass Defect	Relative Mass Defect = $\frac{1e6 \times (exact mass - nominal mass)}{(exact mass)}$
Kendrick Mass Defect	Kendrick Mass Defect = Kendrick mass – nominal Kendrick mass
	where:
	Kendrick Mass = exact mass × nominal mass of Kendrick formula exact mass of Kendrick formula
	Exact mass = Monoisotopic mass of the elemental composition
	Nominal mass = Integer mass
Nominal Mass	Specifies how the node calculates nominal masses.
Rounding	Default: Round
	Selections:
	Floor rounds down.
	Ceiling rounds up.
	 Round rounds to the nearest integer value.
2. Kendrick Formula	as

When you select Kendrick Mass Defect as the Mass Defect Type, this user-specified elemental composition specifies the Kendrick formula.

Formula 1–5	Use to add Kendrick formulas.	
-------------	-------------------------------	--

Compound Class Scoring Node

Use the Compound Class Scoring node to score detected compounds against a set of fragment ions commonly present in the fragmentation scans for a compound class. The node compares the ions (m/z values) detected in the fragmentation scans to the fragments in the selected compound class libraries.

In a processing workflow, connect the Group Compounds node to the Compound Class Scoring node.

The Compound Class Scoring node does the following:

- Annotates the centroids in the fragmentation scans for a compound with the matching fragment structures from the selected compound class libraries.
- Provides a Class Coverage score in the Mass Spectrum view legend.
- Adds the Class Coverage column with the percent coverage to the Compounds table.
- Creates the Compound Class Matches Table—a table related to the Compounds table.

When you add the Compound Class Scoring node to a processing workflow, you must select the compound class fragment lists. For information about adding compound class fragment lists to the Compound Classes library, see "Modifying the Compound Classes List" on page 535.

Table 58 describes the parameters for the Compound Class Scoring node.

Parameter	Description	
1. General Settings		
Compound Classes	Select the compound classes that you want to use for the fragment search.	
	Compound classes contain a list of fragment structures and m/z values that the application compares to the fragmentation scans for each detected compound (see "Modifying the Compound Classes List" on page 535).	
S/N Threshold	Specifies the signal-to-noise threshold for the centroids in the fragmentation spectra. The application ignores centroids below the signal-to-noise threshold. The application attempts to match centroids with m/z values above the threshold to the fragment structures in the selected compound classes.	
	Default: 50	

 Table 58.
 Compound Class Scoring node parameters (Sheet 1 of 2)

Parameter	Description
High Acc. Mass Tolerance	Specifies the mass tolerance for high-resolution mass spectra measured in the Orbitrap mass analyzer of a Thermo Scientific mass spectrometer.
	Default: 2.5 mmu; Minimum: 0.0; Maximum: Unchecked
Low Acc. Mass Tolerance	Specifies the mass tolerance for low-resolution mass spectra measured in the ion trap mass analyzer of a Thermo Scientific mass spectrometer.
	Default: 0.5 Da; Minimum: 0.0; Maximum: Unchecked
Use Full MS Tree	Specifies whether scoring is applied on the full spectrum tree or only the MS2 scans.
Allow DIA Scoring	Specifies whether the node uses DIA scans for scoring when there are no available data-dependent scans. If set to false, the node annotates DIA scans, but it does not use them for scoring.
	Default: True

 Table 58.
 Compound Class Scoring node parameters (Sheet 2 of 2)

Generate Molecular Networks Node (Beta)

Use the Generate Molecular Networks node to determine and visualize the similarity between the various compounds detected by the Detect Compounds node.

Table 59 describes the parameters for the Generate Molecular Networks node.

Table 59. Connect Similar Compounds node parameters (Sheet 1 of 4)

Parameter	Description
1. Spectral Similarity	
These parameters defi similarity score betwee	ne how fragmentation spectra are compared to determine the en two compounds.
Use Full MSn Tree	Specifies whether the node determines the spectral similarity from the full MSn tree or only from the MS2 spectra.
	Default: True
Match Mass Shift	For fragments that are not a direct match by mass, specifies whether the shift to the fragment's mass must match the mass shift between the two compounds.
	Default: True

Parameter	Description	
Match Transformations	For fragments that are not a direct match by mass, specifies whether the shift to the fragment's mass must match the mass of the assigned transformation between the two compounds.	
	Default: True	
Variate Transformations	For fragments that are not a direct match by mass, specifies whether the masses of all the variations of the individual steps of the assigned transformation between the two compounds will be used as the expected fragments shifts. Fragments shifted by these masses will be considered as matching.	
S/N Threshold	Specifies the signal-to-noise threshold for the centroids in the fragmentation spectra. The node ignores centroids below the signal-to-noise threshold—that is, the node attempts to match only centroids with m/z values above the threshold.	
	Default: 50	
Mass Tolerance	Specifies the mass tolerance to use for fragments matching.	
Min. Fragment m/z	Specifies the minimum m/z values of the centroids to consider.	
2. Transformations		
Phase I	Specifies the set of possible single-step Phase 1 transformations.	
	Default: All check boxes are clear.	
Phase II	Specifies the set of possible single-step Phase II transformations.	
	Default: All check boxes are clear.	
Others	Specifies other possible single-step transformations.	
	The node treats Others transformations as Phase I transformations.	
Max. # Phase II	Specifies the maximum number of Phase II steps to be applied.	
	Default: 1; range: 1–10	
Max. # All Steps	Specifies the maximum number of all steps to be applied.	
	All steps that occur as a result of the selections in the Dealkylation area equal one step in the maximum number of all steps—that is after the node applies the steps in the Dealkylation area, the remaining number of possible steps is equal to the Max. # All Steps – 1.	
	Default: 3; range: 1–10	

 Table 59.
 Connect Similar Compounds node parameters (Sheet 2 of 4)

Table 59.	Connect Similar	Compounds node	parameters (Sheet 3 of 4)
-----------	-----------------	----------------	---------------------------

Parameter	Description
-----------	-------------

3. Applied View Filters

These parameters filter out (hide) connections with lower confidence in the Similar Compounds table for a compound. This filtering has no effect on the stored data—that is, you can click the filter icon on the Similar Compounds tab to undo all of these filters.

Require Transformation	When set to True, similar compounds without assigned transformations are hidden.
	Default: True
Require MSn	When set to True, similar compounds without a spectral similarity score are hidden.
Min. MSn Score	Specifies the minimum MSn Score value for a similar compound. If a similar compound has no fragmentation data, the filter is not applied.
Min. MSn Coverage	Specifies the minimum Forward or Reverse Coverage value for a connection. The connection is hidden only if both values are below this threshold. If a connection has no fragmentation data, this filter is not applied.
Min. # Fragments	Specifies the minimum Forward or Reverse Matches value for a connection. The connection is hidden only if both values are below this threshold. If a connection has no fragmentation data, the filter is not applied.

4. Applied Thresholds

These parameters define the filtering rules applied to the stored connections. Compounds that fall below these thresholds are not stored as similar compounds. The main purpose is to filter out low confidence connections to reduce the amount of stored data.

Require Transformation	If set to True, removes connections without assigned transformations.
	Default: True
Require MSn	If set to True, removes connections without a spectral similarity score.
	Default: True
Min MSn Score	Specifies the minimum MSn Score value for a stored connection. If a connection has no fragmentation data, the filter is not applied.

Parameter	Description
Min MSn Coverage	Specifies the minimum Forward or Reverse Coverage value for a stored connection. The connection is removed only if both values are below this threshold. If a connection has no fragmentation data, the filter is not applied.
Min. # Fragments	Specifies the minimum Forward or Reverse Matches value for a connection to be stored (i.e. only if both values are below threshold the connection is removed). If a connection has no fragmentation data, the filter is not applied, and the connection may still be stored.

 Table 59.
 Connect Similar Compounds node parameters (Sheet 4 of 4)

Pattern Scoring Node

Use the Pattern Scoring node to provide a spectrum fit score (SFit%) for each detected compound. The Pattern Scoring node compares the measured isotope pattern for each detected compound to a defined isotope pattern—that is, it compares the mass shifts and intensities of the centroids in the isotope pattern for the detected compound to a set of defined mass shifts and relative intensities. Use the Pattern List Editor dialog box to store defined isotope patterns for your analyses.

The Pattern Scoring node adds the Pattern Matches column to the Compounds table and creates the related Matched Patterns Table.

Table 60 describes the parameters for the Pattern Scoring node.

Parameter	Description	
1. General Settings		
Isotope Patterns	Specifies the isotope patterns to be used for scoring.	
Mass Tolerance	Specifies the mass tolerance for calculated elemental compositions and pattern matching. Default: 5 ppm; range: 0.0–no limit	
Intensity Tolerance [%]	Specifies the relative intensity tolerance of the mass spectral peaks in the isotope pattern. Default: 30; range: 0.01–100.0	
S/N Threshold	Specifies the signal-to-noise threshold for the search. The application ignores isotopes with a theoretical intensity value below this threshold.	

 Table 60.
 Pattern Scoring node parameters (Sheet 1 of 2)

Parameter	Description
Min. Spectral Fit [%]	Specifies the minimum required spectral fit value as a percentage.
	Range: 0 to 100%

Table 60. Pattern Scoring node parameters (Sheet 2 of 2)

Expected Compounds Nodes

These nodes extract information about the compounds that you expect to find in the input file set. The Find Expected Compounds and FISh Scoring nodes require structural information about the targeted compounds. You supply this information by adding it to the Expected Compounds library.

- Find Expected Compounds Node
- FISh Scoring Node
- Generate Expected Compounds Node
- Group Expected Compounds Node
- Mark Background Compounds Node-Expected Compounds
- Merge Features Node

Find Expected Compounds Node

Use the Find Expected Compounds node to search for compounds in the compound ions list provided by one or more Generate Expected Compounds nodes.

Using the input from one or more Generate Expected Compounds nodes, the Find Expected Compounds node looks for expected compounds in the MS1 scans filtered through the data processing nodes. The expected compounds are the parent compounds that the Generate Expected Compounds nodes provide to the Find Expected Compounds node, and the reaction products for these parent compounds. Each Generate Expected Compounds Node predicts the reaction products by using the user-specified Dealkylation step and the user-specified transformation steps. The Dealkylation step can comprise multiple dealkylation and dearylation reactions.

The processing results for the Find Expected Compounds node appear in these tables: Expected Compounds Table, Expected Compounds per File Table, Expected Formulas Table, and Expected Features Table.

For more information about how the application finds expected compounds, see "Targeted Workflows for Expected Compounds" on page 16.

Table 61 describes the parameters for the Find Expected Compounds node.

Parameter	Description
1. General Settings	
Mass Tolerance	Specifies the mass tolerance that the node uses to create each extracted ion chromatogram (XIC).
	Default: 5 ppm; range: 0.1 to 20 ppm
Intensity Tolerance [%]	Specifies the relative intensity tolerance that the node uses for isotope pattern comparison.
	Default: 30%; range: 0–100
Intensity Threshold [%]	Specifies the minimum intensity relative to the base peak for an isotopic peak in an isotope pattern simulation. The application does not search for isotopic peaks below the specified intensity threshold.
	Default: 0.1; range 0.01–10.0
SN Threshold	Specifies the signal-to-noise threshold for the pattern search. Isotopes that are expected to be below this intensity level are not required. This means that if the intensity for an isotope in the theoretical isotope pattern (based on the expected elemental composition) is below the signal-to-noise threshold, the application ignores the isotope and requires one less matching isotope in the measured isotope pattern.
	Range: 1 or greater
Min. # Isotopes	Specifies the minimum number of isotopes (mass spectrum peaks in a centroided mass spectrum) that must match the theoretical isotope pattern of the expected elemental composition.
	Default: 2; range: 1 to no limit
Minimum Peak Intensity	Specifies the minimum apex intensity, in counts, of the detected chromatographic peak. The node discards chromatographic peaks below this intensity threshold.
	Default: 1000; Minimum: 0.0

 Table 61.
 Find Expected Compounds node parameters (Sheet 1 of 2)

Parameter	Description
Average Peak Width	Specifies the average chromatographic peak width (FWHM) in the filtered time range.
	Default: 0 (automatic peak width detection); range: unchecked
	When this value is set to 0, the node automatically determines the average peak width.
	IMPORTANT The node detects no chromatographic peaks in the following cases:
	• The filtered retention time is too small compared to the determined or user-specified average peak width value. For information about filtering the scan data, see "Select Spectra Node" on page 174.
	• The determined or user-specified average peak width value is too small compared to the scan rate of the instrument.
	For example, if the instrument acquires a full (MS1) scan every 0.01 minutes, do not enter an Average Peak Width value of less than 0.02 (2×0.01 minutes), as the peak detection algorithm requires a minimum of three data points to detect a chromatographic peak.
	For best results, keep the default setting of 0, which turns on automatic peak width detection. Enter a nonzero value only when the automatically detected peak width is not suitable or fails for your chromatographic method.

Table 61. Find Expected Compounds node parameters (Sheet 2 of 2)

FISh Scoring Node

Use the FISh Scoring node to provide a confirmation score for compounds that the Find Expected Compounds node detects and to annotate the fragmentation spectra for these compounds. The FISh Scoring node requires data-dependent fragmentation (DDF) scans to calculate the FISh coverage scores for related structures.

For information about how the node calculates the confirmation score, see "FISh Scoring for Targeted Compounds and Proposed Structures" on page 36.

Table 62 describes the parameters for the FISh Scoring node.

Table 62. FIS	h Scoring node	parameters	(Sheet 1 of 2)
---------------	----------------	------------	----------------

Parameter	Description
1. General Settings	
Annotate Full Tree	Specifies whether the node annotates the full spectrum tree or only the MS2 scans in the Mass Spectrum view.
	For information about viewing the FISh annotations in the Mass Spectrum view, see "Working with the Mass Spectrum View" on page 305.
	Default: True
Match Transformations	Specifies whether the node matches fragments with transformation shifts.
	Default: True
S/N Threshold	Specifies the signal-to-noise threshold for centroids. The node ignores centroids below this threshold in the fragmentation (MS/MS, MS ³ , and so on) spectra.
	Default: 3
High Acc. Mass Tolerance	Specifies the mass tolerance for high-resolution mass spectra measured in the Orbitrap mass analyzer of a Thermo Scientific mass spectrometer.
	Default: 2.5 mmu; Minimum: 0.0; Maximum: Unchecked
Low Acc. Mass Tolerance	Specifies the mass tolerance for low-resolution mass spectra measured in the ion trap mass analyzer of a Thermo Scientific mass spectrometer.
	Default: 0.5 Da; Minimum: 0.0; Maximum: Unchecked
2. Fragment Prediction Se	ttings
Use General Rules	Specifies whether the node uses the general fragmentation rules.
	Default: True
Use Libraries	Specifies whether the node uses fragmentation libraries for fragment prediction.
	Default: True
	Note Using fragmentation libraries to predict fragments adds significant time to data processing; however, it also provides significantly more predicted fragments.

Parameter	Description
Max. Depth	Specifies the maximum number of steps in the fragmentation pathway.
	Range: 1 to 20
Aromatic Cleavage	Specifies whether the node includes a cleavage step in the fragmentation pathway for highly aromatic structures—that is, for aromatic structures where n in Huckel's rule is 2 or higher.
	4n + 2 = 0 or a positive integer for the number of electrons in the delocalized, conjugated p-orbital cloud
	For example, the following structure is aromatic with an n value of 2.
	HONN
	Default: True
Min. Fragment <i>m/z</i>	Specifies the minimum <i>m/z</i> value of a fragment ion to be generated by the prediction fragmentation pathway.
	Default: 50

Table 62. FISh Scoring node parameters (Sheet 2 of 2)

Generate Expected Compounds Node

Use the Generate Expected Compounds node to generate a list of m/z values for the ionized compounds that you expect to find in a sample. The list includes the parent compounds and their possible dealkylation, dearylation, and transformation products. The application generates the list by using the structures of the parent compounds, the user-specified transformation lists and number of combinatory steps, and the user-specified ionic species.

The default transformations library contains common Phase 1 and Phase 2 metabolic transformations. If the transformation list does not include the possible transformations for your compound, add them to the transformations library as described in "Adding or Editing Transformations with the Transformation Editor" on page 512.

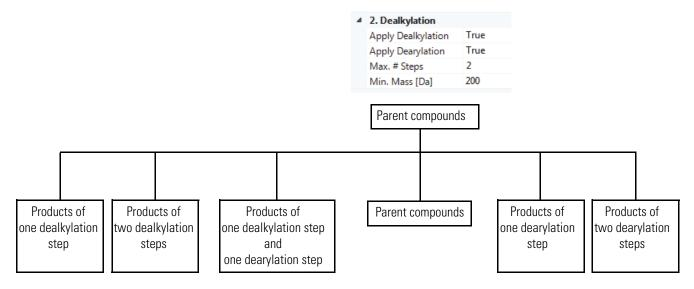
When you add transformations to the library, you can assign them to one of the following groups: Phase I, Phase II, or Others. The Generate Expected Compounds node treats transformations assigned to the Others group as Phase I transformations; that is, it applies Phase I and Others transformations before it applies Phase II transformations.

To predict the transformation products for the selected parent compounds, the Generate Expected Compounds node follow these rules:

1. When the user enables Dealkylation, apply the dealkylation steps first. If a subsequent transformation reverses the dealkylation step, reject the subsequent transformation. When the user enables both Dealkylation and Dearylation, apply both of these steps first, and then determine two separate reaction pathways for the remaining transformation steps.

Consider all steps under Dealkylation together as one step. For example, consider the selections shown in Figure 70 as one step in the total set of reaction pathways and create separate reaction pathways. Apply the transformation steps on the parent compound and the reaction products from the dealkylation pathways.





- 2. When more than one reaction pathway produces the same elemental composition, use the pathway with the lowest number of transformation steps.
- 3. Reject transformations that remove elements that are not present. For example, do not apply an oxidative dechlorination step if the compound does not contain chlorine.
- 4. For Phase I and Others transformations, limit the maximum number of times to apply the transformation on a single compound to the lower of these two values:
 - Max Occurrence setting for the transformation in the Transformations library

For example, for the oxidation transformation, the default value for Max Occurrence is 3.

Name		Arrivin	g Group	Arriving	g Modification	ΔΜ [Da]	Phase	Max Oc	currence 🔻
Aa	•	<u>A</u> a	*	<u>A</u> a	*	=	•	<u>A</u> a	=	•
Oxidation		0		0		15.	99491	Phase1		3

 Maximum number of steps specified by the node's Max. # All Steps parameter minus any previously applied Dealkylation step.

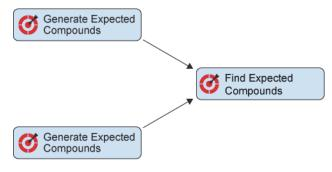
- 5. For Phase II transformations, limit the maximum number of times to apply the transformation on a single compound to the lowest of these three values:
 - Max Occurrence setting for the transformation in the Transformations library
 - Maximum number of steps for all reactions (setting for node's Max. # All Steps parameter) minus any previously applied Dealkylation or Transformation step
 - Maximum number of steps for a Phase II transformation (setting for node's Max. # Phase II parameter)

You can connect one or more Generate Expected Compounds nodes to the Find Expected Compounds Node and the Filter By Mass Defect Node.

Table 63 describes the parameters for the Generate Expected Compounds node. The application cannot use the processing workflow until you select a compound from the Compound list.

If you want to generate expected compounds for more than one parent compound, do the following:

• To target multiple compounds with a different set of transformation rules for each compound, add multiple Generate Expected Compounds nodes to the processing workflow.



• To target multiple compounds with the same set of transformation rules, add one Generate Expected Compounds node to the processing workflow and select multiple compounds in the node's Compound list.

Par	ameters of 'Generate Expe	ted Compounds'	
Sh	ow Advanced Parameters		
⊿	1. Compound Selection		
	Compound	Caffeine (C8 H10 N4 O2); Omeprazole (C17 H19 N3 O3 S)	-
4	2. Dealkylation Apply Dealkylation Apply Dearylation	Show Checked Only (2/3) Filter	
	Max. # Steps Min. Mass [Da]	 2-Hydroxy-4-quinolincarboxylic acid (C10 H7 N O3) Caffeine (C8 H10 N4 O2) 	*
4	3. Transformations Phase I	Omeprazole (C17 H19 N3 O3 S)	-
	Phase II Others	Check All Uncheck All	
	Max. # Phase II Max. # All Steps		
⊿	4. Ionization		
	Ions	[M+H]+1	

Table 63. Generate Expected Compounds node parameters (Sheet 1 of 3)

Parameter	Description	
1. Compound Selection		
Compound	Specifies the parent compounds that the node uses to build a list of possible product compounds.	
Show Checked Items Only	Displays only the selected compounds.	
Check All	Selects all compounds in the list.	
Uncheck All	Undoes all selections.	
transformation produ Expected Compound 2. Dealkylation	cts of specific compounds, you must first add the compounds to the s library.	
Apply Dealkylation	When you select True, the node applies the dealkylation transformations for the specified compound before applying other transformations. Default: True	
Apply Dearylation	When you select True, the node applies the dearylation transformations for the specified compound before applying other transformations.	
	Default: False	

Parameter	Description			
Max. # Steps	Specifies the maximum number of Dealkylation steps.			
	For example, if you select True for Dealkylation, True for Dearylation, and 1 for the Max. # Steps, the node applies up to one dealkylation step and up to one dearylation step as the initial Dealkylation step in the set of reaction pathways. For another example, see Figure 70 on page 235.			
	Parent compound			
	Products of one dearylation stepProducts of one dearylation step			
	Default: 1; Selection: 1–10			
Min. Mass [Da]	Specifies the minimum mass of the dealkylation product.			
	Default: 200			
3. Transformations				
Phase I	Specifies the set of possible Phase 1 transformations.			
	Default: All check boxes are clear.			
Phase II	Specifies the set of possible Phase II transformations.			
	Default: All check boxes are clear.			
Others	Specifies other possible transformations.			
	The node treats Others transformations as Phase I transformations.			
Max. # Phase II	Specifies the maximum number of Phase II steps to be applied.			
	Default: 1; range: 1–10			

Table 63. Generate Expected Compounds node parameters (Sheet 2 of 3)

Parameter	Description
Max. # All Steps	Specifies the maximum number of all steps to be applied.
	All steps that occur as a result of the selections in the Dealkylation area equal one step in the maximum number of all steps—that is, after the node applies the steps in the Dealkylation area, the remaining number of possible steps is equal to the Max. # All Steps – 1.
	Default: 3; range: 1–10
4. Ionization	
Ions	Specifies the possible ionic species.
	Default: [M+H]+1 (protonated species for the positive mode)
	Note The ion definitions library that the application provides contains the common ionic species associated with the positive and negative modes for the electrospray ionization-mass spectrometry (ESI-MS) technique. If the Ions list does not include the possible ionic species for your analysis, add the ion definition to the Ion Definition library as described in "Modifying the Ion Definitions List" on page 505.

 Table 63.
 Generate Expected Compounds node parameters (Sheet 3 of 3)

Group Expected Compounds Node

Use the Group Expected Compounds node to combine similar components (chromatographic peaks with the same MW×RT dimensions) that the Find Expected Compounds node finds across the input file set. This node combines chromatographic peaks by using their chemical formula (resulting from the dealkylations and dearylations and transformations of a parent compound) and retention time. This node also selects the best representative MS1 scan and fragmentation tree, which other nodes, such as Predict Compositions, Search mzCloud, Search mzVault, and so on, use for identification.

Table 64 describes the parameters for the Group Expected Compounds node.

Table 64.	Group Expected	Compounds node	parameters (Sheet 1 of 2)
-----------	----------------	----------------	---------------------------

Parameter	Description	
1. General Settings		
RT Tolerance [min]	Specifies the retention time tolerance, in minutes, that the node uses to group mass peaks generated from the same parent compound through the same reaction pathway.	
	Default: 0.05; range: 1.0	

Parameter	Description
2. Fragment Data Sel	ection
Preferred Ions	Select the preferred ions (adducts) from the list. The application uses the list to select the best fragmentation data for each compound to submit to an mzCloud or mzVault search.

 Table 64.
 Group Expected Compounds node parameters (Sheet 2 of 2)

Mark Background Compounds Node–Expected Compounds

Use the Mark Background Compounds node to flag compounds that are also found in the sample blanks (Sample Type—Set to Blank). The Mark Background Compounds node for a targeted analysis has a blue background.



For information about editing the assigned sample types, see "Editing the Sample Type and Study Factor Values" on page 110.

Table 65 describes the parameters for the Mark Background Compounds node.

Table 65. Mark Background Compounds node parameters (Sheet 1 of 2)

Parameter	Description
1. General Settin	gs
Max. Sample/Blank	Below this ratio threshold, the node labels the expected compound as a background compound in the Expected Compounds table. If the input file set includes more that one blank sample, the node uses the largest peak area for the compound in the blank input files as the peak area of the blank. $\frac{\text{Peak Area}_{\text{Sample}}}{\text{Peak Area}_{\text{Blank}}} < \text{Specified value}$
	Default: 5
	When this setting is 0, the node ignores this parameter.

Parameter	Description
Max. Blank/Sample	Below this ratio threshold, the node labels the expected compound as a background compound in the Expected Compounds table. If the input file set includes more that one blank sample, the node uses the largest peak area for the compound in the blank input files as the peak area of the blank. $\frac{\text{Peak Area}_{\text{Blank}}}{\text{Peak Area}_{\text{Sample}}} < \text{Specified value}$
	If the compound is found in a blank sample but not in a non-blank sample, (ratio X/0), the node marks the compound as a background compound.
	Default: 0 (The node does not use this parameter to mark background compounds.)
Hide Background	When the Hide Background parameter is set to True, the tab for the Expected Compounds Table includes a filter icon (Expected Compounds radius and the expected compounds that fall below the threshold are hidden. Clicking the filter icon displays the filtered compounds.
	When the Hide Background parameter is set to False, the background compounds appear in the result table.
	The Background column is a hidden column in the Expected Compounds table. When a compound is flagged as a background compound, its Background check box is selected.
	Background
	Default: True

Table 65. Mark Background Compounds node parameters (Sheet 2 of 2)

Merge Features Node

Use the Merge Features node to do the following:

- Combine the expected compounds found by the Find Expected Compounds node and the unknown compounds found by the Detect Compounds node by using their chromatographic retention time and *m/z* values.
- Create the Merged Features table that includes four status columns: Ion Conflict Status, Detect Compounds, Find Expected Compounds, and Custom Explanations.
- Link the manual peaks table to the related Compounds table.

Table 66 describes the parameters for the Merge Features node.

Parameter	Description	
1. General Settings		
Mass Tolerance	Specifies the mass tolerance for chromatographic peak grouping.	
	Default: 5 ppm; range: 0.1–20 ppm	
RT Tolerance [min]	Specifies the retention time tolerance, in minutes, for chromatographic peak grouping.	
	Default: 0.05; range: 0.0–1.0	

Table 66. Merge Features node parameters

Post-Processing Nodes

These nodes provide additional information about the detected compounds.

- Descriptive Statistics Node
- Differential Analysis Node
- Export Xcalibur Inclusion or Exclusion List Node
- Scripting Node

Descriptive Statistics Node

Use the Descriptive Statistics node to store descriptive statistics for each expected or unknown compound in the result file and add the following hidden columns to the Compounds and Expected Compounds tables: Mean Area, Median Area, Minimum Area, Q1 Area, and Q3 Area.

To store the descriptive statistics in the result file, the processing workflow must include the Group Expected Compounds node for targeted workflows and the Group Compounds node for untargeted workflows.

The Descriptive Statistics node has no parameters.

Note The Descriptive Statistics node generates the descriptive statistics for individual compounds; it has no effect on the Descriptive Statistics view that is available for a result file.

Differential Analysis Node

Use the Differential Analysis node to calculate the statistics for a differential analysis (fold change, ratio, p-values, and so on), store this data in the result file, and create a volcano plot in the Differential Analysis view by using the data stored in the Compounds Table or Expected Compounds Table. A volcano plot is a type of scatter plot for replicate data where the *x* axis represents the log₂ of the fold change between two sample groups (generated ratio), and the *y* axis represents the negative log₁₀ of the p-value (test of significance) of the fold change.

This node requires input from the Group Compounds node or the Group Expected Compounds node for a sample set with replicate data points and generated ratios. If the Grouping and Ratios page of an analysis does not contain generated ratios, the following confirmation message appears:

No quan ratios defined in 'Grouping & Ratios' tab

Table 67 describes the parameter for the Differential Analysis node.

Parameter	Description			
1. General Settings				
Log10 Transform Values	Specifies whether to calculate \log_{10} values of the chromatographic peak areas before storing the data in the result file.			
	Default: True			

Export Xcalibur Inclusion or Exclusion List Node

Use the Export Xcalibur Inclusion/Exclusion List node to create an Inclusion/Exclusion mass list for a Thermo Scientific mass spectrometer.

Table 68 describes the parameters for the Export Xcalibur Inclusion or Exclusion List node.

Parameter	Description			
1. General Settings				
File Name	Specifies the file name for the inclusion/exclusion list.			
Selected Instrument	Bases the format of the inclusion/exclusion list on the selected MS.			
	Selections: LTQ Orbitrap, Orbitrap Fusion, Q Exactive			
2. Use Filter Set				
Add Filter Set	Specifies the filters sets (FILTERSET file type) that you want to apply to the generated inclusion list or exclusion list.			

Table 68. Export to Xcalibur Inclusion/Exclusion List node parameters (Sheet 1 of 2)

Parameter	Description				
3. Advanced Settings					
Left RT [min]	Specifies the window to the left of the specified retention time for a mass.				
	Default: 1 min; range: 0.001 to 1000 min				
Left RT [min]	Specifies the window to the right of the specified retention time for a mass.				
	Default: 1 min; range: 0.001 to 1000 min				
Include Isotopic Peaks	Specifies whether to include isotopic peaks in the list.				
	Default: False				
4. LTQ Orbitrap Settings					
Maximum Concurrent Entries	Specifies the maximum number of entries with overlapping time windows.				
	Default: 500; range: 1 to 2000				
Mass Precision Decimals	Specifies the required number of decimal places for the mass values.				
	Default: 5				

 Table 68.
 Export to Xcalibur Inclusion/Exclusion List node parameters (Sheet 2 of 2)

Scripting Node

Use the Scripting Node to perform custom post-processing actions on the data in the result tables.

Tip For information about using the Scripting Node, go to the Resources page of the following web site: https://mycompounddiscoverer.com/.

Table 68 describes the parameters for the Scripting Node.

Table 69. Scripting Node parameters (Sheet 1 of 2)

Parameter	Description		
Executable and Parameters			
Path to Executable	Specifies the path to the executable file, for example, c:\Python37\python.exe.		
Command Line Arguments	Specifies command line arguments provided to the executable, for example, "d:\My Scripts\script.py \" %NODEARGS%.		

Parameter	Description				
Requested Tables and Columns	Specifies the requested result table columns for the executable.				
	If you specify a table without specifying specific columns in the table, the node exports the entire table.				
	Use a colon to separate the table name from the table columns, a comma to separate the table columns, and a semicolon to separate tables.				
	Format:				
	TableName1:Column1, Column2;TableName2:Column3, Column4				
	✤ To enter this list				
	1. Click the browse icon.				
	The Edit Parameter Text for Requested Tables and Columns dialog box opens.				
	2. Do one of the following:				
	• Enter the tables and columns requested. Use a new line for each table.				
	• Click Load File , select a text file that contains the table and column information, and click Open .				
	3. Click OK .				
Use R-Friendly	Specifies whether the column names are R-Friendly.				
Columns	Default: True				
Archive Datafiles	Specifies whether the node creates an archive containing the JSON files and the text (.txt) files.				
	Default: False				

Table 69. Scripting Node parameters (Sheet 2 of 2)

7 Workflow Nodes Post-Processing Nodes

8

Reviewing the Analysis Results

The Compound Discoverer application stores the results of an analysis in a result file (cdResult).

These topics describe how to open result files, filter the result tables to remove non-pertinent data, edit the compound annotations, propose custom structures, and modify and store the layout of the current result file. They also describe how to export mass lists and spectral data.

Contents

- Opening, Closing, and Updating Result Files
- Factory Default Layout for a Result Page
- Modifying the Result Page Layout
- Saving, Restoring, and Managing Layouts
- Editing Compound Annotations
- Adding and Deleting Proposed Structures for a Compound
- Replacing an Annotation with a Structure Proposal
- Applying FISh Scoring
- Using Result Filters for Data Reduction and Creating Filter Sets
- Viewing the Result Summary
- Exporting the Tabular Results to an External File
- Exporting Spectral Data to a New or Existing mzVault Library
- Exporting Compounds to a New or Existing Mass List
- Copying or Saving Graphical Views for Publication

Related Topics

- Chapter 9, "Working with the Graphical Result Views."
- Chapter 10, "Reviewing the Tabular Result Data."

Opening, Closing, and Updating Result Files

During an analysis, the Compound Discoverer application processes a set of input files (Xcalibur RAW files) by using a processing workflow and stores the processing results in a result file (CDRESULT).

When you open a result file for the first time, you see a tabbed document with the default layout in the application window. You can modify the layout and save these changes with the result file. The next time you open the result file, it will open with your custom layout.

To open a result file from a previous version of the application, you must update the file to the current version.

For instructions on how to open, close, and update result files, see these topics:

- Opening Result Files Created in the Current Application Version
- Opening Result Files Created in Previous Versions of the Application
- Update Modes for Legacy Result Files
- Closing a Result File

Opening Result Files Created in the Current Application Version

You can open a result file from the application window, the Start Page, the Job Queue page, or the Analysis Results page. A result file opens as a tabbed document. The tab displays the file name of the result file (*File Name* X).

* To open a result file

- From the application window, do one of the following:
 - From the menu bar, choose File > Open Result. In the Open dialog box, browse to the appropriate folder, select the result file of interest, and click Open.
 - From the menu bar, choose File > Recent Results > recent result file.
- From the Start Page, do one of the following:
 - Under Recent Results, click the blue hyperlink for the result file of interest.
 - Under What Would You Like to Do?, click **Open Result**.
- From the Job Queue page, do one of the following:
 - Double-click the table row for a completed job.
 - Select the table row of a completed job and click **Open Results**.
- From the Analysis Results page, do one of the following:
 - Double-click the table row for a completed analysis.
 - Select the file of interest and click **Open Results**.

Tip You can also drag and drop result files (CDRESULT) from Windows Explorer into the application window.

You can open as many result files as you want. To view a particular result file, click its tab.

Opening Result Files Created in Previous Versions of the Application

* To open a result file from a previous version of the application

1. From the menu bar, choose **File > Open Result**. In the Open dialog box, browse to the appropriate folder, select the result file of interest, and click **Open**.

The File Update Required dialog box opens (Figure 71).

Figure 71. File Update Required dialog box

File Update Required			
The file 'E&L_Oring.cdResult' needs to be updated. How do you want to perform the update?			
🖉 Keep Original File			
☑ Failsafe Update			
Update Cancel			

- 2. Select one of the following update processes:
 - To retain the original result file, select the Keep Original File check box.

The application automatically selects the Failsafe Update check box as well.

• To automatically delete the original file after the update process ends, select only the **Failsafe Update** check box.

When the Failsafe Update check box is selected, the application automatically backs out of the update process if an error occurs.

-or-

• To minimize the processing time, clear both check boxes.

The application immediately processes the file without first creating a temporary file. Errors that occur during the update process can corrupt the file.

IMPORTANT To prevent loss of data when not using the fail-safe mode to update legacy result files, Thermo Fisher Scientific recommends that you manually store copies of these files in another directory.

3. To start the update process, click Update.

Update Modes for Legacy Result Files

When you attempt to open a legacy result file, the application prompts you to select an update process:

- Selecting the Keep Original File check box runs the update process in the fail-safe mode and renames the original result file with an appended version number. This option takes the most processing time, but it prevents data corruption of the original file.
- Selecting the Failsafe Update check box runs the update process in the fail-safe mode, but it does not save the original result file. If the update process fails, the application retains the original result file. You can make another attempt to update the file or you can open the file in a previous version of the application.

Note You can install multiple versions of the Compound Discoverer application on the same data system computer.

• Clearing both check boxes turns off the fail-safe mode. This option takes the least amount of processing time, but it risks the possibility of corrupting the original file and making it unrecoverable.

In the fail-safe mode, the application does the following:

- 1. Creates a temporary file.
- 2. Runs the updates on the temporary file.
- 3. After completing the update process successfully, it does the following:
 - a. Appends the application version to the file name of the original file. Because the application does not reprocess the original file, the file retains its original time stamp.
 - b. Changes the file name extension of the temporary file to CDRESULT. The time stamp for the updated file corresponds to the completion of the update process.

If the update process fails, the application does not rename the original file.

Closing a Result File

You can close a result file (CDRESULT) in two ways.

To close a result file

• Right-click the tab and choose Close.

-or-

• Click the close icon on the document's tab (\boxtimes) .

If the tab is not visible, click the Current Tabs icon, , and select the result file from the list (see "Working with the Tabbed Documents" on page 41).

Factory Default Layout for a Result Page

When you open a result file (see Opening, Closing, and Updating Result Files), it appears as a tabbed page in the application window.

The factory default layout for a result page includes the following items:

- A tab with the result file name
- A Chromatograms view on the top left that is populated with XIC traces for the component with the largest chromatographic peak area across the input files—that is, the view is populated with the XIC traces for the compound listed in the first row of the Compounds table or the Expected Compounds table. The view is zoomed to the start and end points of the chromatographic peak.
- A Mass Spectrum view on the top right that is populated with the MS1 scan with the highest resolution and highest intensity related to the preferred ions across the input files. The spectrum tree to the left includes the MS1 scans and the available fragmentation scans within the following time range for a compound:
 - Peak apex (RT) ± the peak's full width at half maximum (FWHM)

-or-

- Start and end points of the chromatographic peak, as determined by the peak detection algorithm
- A set of tabbed main tables below the two graphical views
- A collapsed area for the related tables below the main tables

Figure 72 shows a result page for a result file generated by using the MetID w Stats Expected w Background Removal workflow template.

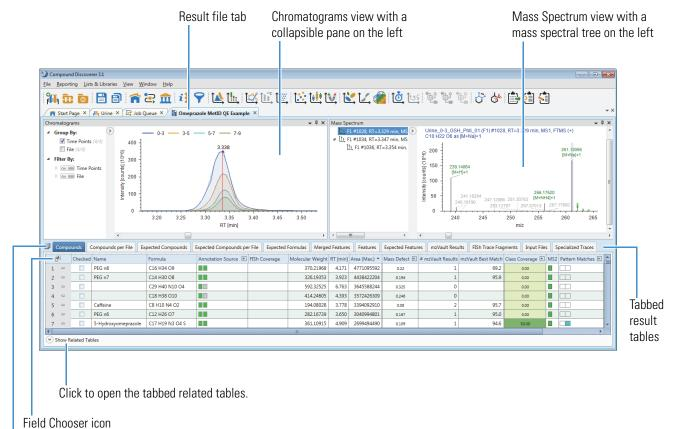


Figure 72. Default layout for a result page

Select Table Visibility icon

You can change the layout as follows:

- The views that you want to display and their location
- The main and related tables and table columns that you want to display and the order of the table columns from left to right
- The data (table rows) displayed or hidden in the result tables

Related Topics

• Opening, Closing, and Updating Result Files

Modifying the Result Page Layout

The result file layout includes the relative positions of the graphical views, the visible result tables, the column arrangement and column fixing if it is enabled, and the applied result filters.

These topics describe some of the ways that you can change the layout of a tabbed page for a result file:

- Floating a Result Page View
- Enlarging a Result Page View
- Collapsible Pane Options for Filtering, Grouping, Coloring, and Discriminating By
- Using a View's Collapsible Pane to Filter the Data By the Study Variable Values
- Using a View's Collapsible Pane to Group the Data
- Showing or Hiding Result Tables

Floating a Result Page View

* To change a docked view on a result page to a floating view

Do one of the following:

- Right-click the view's title bar and choose Floating.
- Double-click the view's title bar.
- -or-
 - Drag the view away from its docked position.

Enlarging a Result Page View

* To enlarge a floating view for a result page to fill the screen

Double-click the view's title bar.

To restore the previous size of a floating window, double-click the view's title bar again.

Collapsible Pane Options for Filtering, Grouping, Coloring, and Discriminating By

These four views on a result page include a collapsible pane of filtering options and grouping, coloring, or discriminating by options:

- Chromatograms view (see "Working with the Chromatograms View" on page 294)
- Trend Chart view (see "Working with the Trend Chart View" on page 332)

- Principal Component Analysis view (see "Working with the Principal Component Analysis View" on page 341)
- Partial Least Squares Discriminant Analysis (see "Working with the Partial Least Squares Discriminant Analysis View" on page 358)
- Descriptive Statistics view (see "Working with the Descriptive Statistics View" on page 347)

Note If the analysis does not include samples with different study factor values or sample types, the application cannot group the samples, and only the Files check box appears in the Group By list.

Table 70 describes the effect of clearing and selecting the check boxes in the collapsible pane at the left of a view.

Objective	Selection
Remove or display data points	Use the check boxes under Filter By to remove or display data points by sample type, study factor value, or individual file. By default, the application selects all the check boxes with the exception of Blank sample types
Visually distinguish data points by grouping them	For the Chromatograms view, use the check boxes under Group By to colorize the traces by group. The application duplicates the check boxes in the Study Variables area of the Grouping and Ratios page of the analysis. For the Trend Chart view, use the check boxes under Group By to change how the data points for the selected compound are grouped across the input files.
Distinguishing data point by color	For the Principal Component Analysis view, use the Color By check boxes to distinguish the principal components by color. For the Descriptive Statistics view, use the Color By check boxes to visually group the box plots by color.
Discriminate by specific study variables	For the Principal Least Squares–Discriminant Analysis view, use the Discriminate By check boxes to select the study variables for the supervised analysis.

Table 70. Effect of clearing or selecting the check boxes in the collapsible pane

Table 71 describes the Group By and Filter By options in the collapsible pane.

Feature	Description			
(<i>X</i> / <i>Y</i>)	The left integer is the number of samples that are selected under Filter By (and that also contain the selected compound). The right integer is the number of samples that contain the selected compound.			
ON/OFF toggle for	ON—The check boxes are available.			
the Filter By items	OFF—The check boxes are unavailable and the items are not filtered out.			
Check boxes				
Group By <i>Study</i> Variable	Selecting one or more of these check boxes groups the samples with the same value or values for the selected study variable or variables and displays the groups in different colors.			
	By default, the application duplicates the selection in the Study Variables pane on the Grouping and Ratios page of the analysis. Study variables include the study factor values and the sample types.			
Group By Sample	Selecting this check box displays the data points for the selected samples in the Filter By area in different colors.			
	Default: Clear			
Filter By Study Factor	Select these check boxes to display data for one or more study factors.			
	By default, the application selects all the study factors.			
Filter By Sample Type	Select these check boxes to display data for one or more sample types.			
	By default, for the Chromatograms view, the application selects all sample types. For the statistical views, the application excludes the Identification Only sample type and clears the Blank sample type check box.			
Filter By File	Select these check boxes to display data from one or more of the files.			
	By default, the application selects all of the files.			

Table 71. Options in the collapsible left pane

Using a View's Collapsible Pane to Filter the Data By the Study Variable Values

You can filter the data points displayed in a result page view by selecting or clearing the check boxes for the study variable values in the collapsible pane.

* To filter data points by the study variable values

- 1. If the view's collapsible pane is closed, click the icon, ((), in the upper left corner of the view.
- 1. Under Filter By, click the expand icon to the left of the study variable name to open the values list.
- 2. Clear the check boxes for the items that you want to hide or values that you want to remove from the statistical calculations.

Using a View's Collapsible Pane to Filter the Data By Input File

You can filter the data points displayed in a result page view by selecting or clearing the check boxes for the input files in the collapsible pane.

To filter the data by selected files

- 1. Under Filter By, click the expand icon to the left of File to open the File list.
- 2. Clear the check boxes for the files that you want to exclude from the display or the statistical calculations.

Note By default, for the statistical views, the check box for the Blank sample type under Filter By is clear.

Using a View's Collapsible Pane to Group the Data

You can group the data points displayed in a result page view by selecting or clearing the check boxes under Group By in the collapsible pane.

To group the data points in a view

- 1. To group the data by the study variables or by the individual files, select a row in the active result table. Then, select one or more check boxes under Group By
- 2. To view a color legend of the sample groups, right-click the Chromatograms view and choose **Display Options > Show Legend**.

Showing or Hiding Result Tables

The set of result tables in a result file depends on the processing workflow.

By default, some of the result tables are hidden. For example, the Adducts table is hidden for both targeted and untargeted analyses and the Filled Gaps related table is hidden for an untargeted analysis.

✤ To show or hide result tables

- 1. Open a result file (see "Opening, Closing, and Updating Result Files" on page 248).
- 2. Click the **Select Table Visibility** icon, I, to the left of the result table tabs.

The Select Visible Tables dialog box opens.

3. Select the check box that corresponds to the table that you want to show, or clear the check box that corresponds to the table that you want to hide.

By default, the Adducts and Filled Gaps tables are hidden.

4. Click **OK** to accept the changes.

Saving, Restoring, and Managing Layouts

The layout of the result file includes the location of the graphical views, the result tables that you want to display, the columns and rows that you want to display in the result tables, the filter set, the custom annotations, and the Group By and Filter By settings.

In addition to the factory default layout, the application comes with the following custom layouts:

- Identification—Adds the Structure column to the Compounds and Expected Compounds tables. Opens the primary set of related tables to the Structure Proposals table. Closes the search result tables if they are visible.
- Quantification—Opens the Trend Chart view. Opens the primary set of related tables to the Structure Proposals table. Closes the search result tables if they are visible.
- Stable Isotope Labeling—Opens the Isotopologues Distribution Chart, the Trend Chart, and the Metabolika Pathways view. In the Compounds table, hides the following columns: #Metabolika Pathways, Avg. Exchange, FISh Coverage, and Metabolika Pathways,
- Statistics—Opens the Differential Analysis, Trend Chart, Principal Component Analysis, and Partial Least Squares Discriminant Analysis views. Closes the Chromatograms and Mass Spectrum views. Closes the search result tables if they are visible.

Follow these topics as needed:

• Saving the Current Layout of a Result File

- Resetting the Layout to the Factory Defaults
- Creating a Custom Layout
- Applying a Custom Layout
- Managing the Layouts

Saving the Current Layout of a Result File

* To save the current layout of the result file

With the result file selected as the active page, do one of the following:

• In the toolbar, click the Save the Currently Active Item icon, 📙.

-or-

• From the menu bar, choose **File > Save** (save the currently active item).

Resetting the Layout to the Factory Defaults

* To reset the layout to the factory default settings

With the result file selected as the active page, from the menu bar, choose **Window** > **Reset Layout**.

The application closes the result file, and then reopens the result file to the factory default layout.

Creating a Custom Layout

- ✤ To create a custom layout
- 1. Open a result file and modify its layout.

For details, see "Opening, Closing, and Updating Result Files" on page 248 and "Modifying the Result Page Layout" on page 253.

2. From the menu bar, choose **Window > Save Layout**.

The Save Result Layout dialog box opens.

Save Result Layout	
Layout Name:	
New Layout 1	
	<u>O</u> K <u>Cancel</u>

3. Name the layout and click **OK**.

Applying a Custom Layout

✤ To apply a custom layout

- With the result file selected as the active page, from the menu bar, choose Window > Apply Layout.
- 2. Select a layout from the list or use the hot keys.

Managing the Layouts

✤ To manage the layouts

1. From the menu bar, choose **Window > Manage Layouts**.

The Manage Result File Layouts dialog box opens. The list order corresponds to the hot keys in the Apply Layout list. For example, for this list, the hot key combination for the Statistics layout is CTRL+ALT+1.

Manage Result File Layouts	×
Saved Layouts: Identification Quantification Stable Isotope Labeling Statistics Rename Delete As Default Close	Use these buttons to change the order of the layouts. Changing the order changes the associated hot keys. The first 10 layouts in the list have associated hot keys.

- 2. Do the following as needed:
 - To delete a layout, select it and click **Delete**.
 - To rename a layout, select it and click **Rename**. Then, in the Rename Result Layout dialog box, rename the layout and click **OK**.
 - To make the layout the default layout, select it and click As Default.
 - To change the list order, select a layout, and use the Up/Down buttons, , to move the layout up or down in the list.

Editing Compound Annotations

Annotations include the compound name, formula, annotation source, FISh coverage score, and structure.

Use the Compound Annotation Editor dialog box to edit the annotations for compounds of interest in the Compounds Table, Expected Compounds Table, and the Structure Proposals Table. Saving a custom annotation overwrites the original processing results.

✤ To edit a compound annotation

1. Double-click the row of interest in the Compounds table, Expected Compounds table, or related Structure Proposals table.

The Compound Annotation Editor dialog box opens (Figure 73).

The application automatically populates the Molecular Weight/Error in Da, and Molecular Weight to Fit boxes. If the formula, structure, and name are available, the application also populates these fields.

- 2. To add a structure to the drawing area, do any of the following:
 - Use the structure drawing tools. See "Using the Structure Drawing Tools or Commands" on page 542.
 - Open a structure file. See "To open a structure file" on page 501.
 - Run a ChemSpider search and select one of the hits. See "Finding a Structure in the ChemSpider Database" on page 541.

3. Click **Save** to save your custom annotations in the result file.

Figure 73. Compound Annotation Editor dialog box

Open (structure file) Save	Compound Annotation	Editor $\checkmark \land \# \# \downarrow^r \textcircled{O}$ H H H H H H H H		Drawing area
(structure file)				
	Description FISh Scoring			
The application	Formula:		Formula to fit:	
	C13 H14 N2 O2			
	Molecular weight / error in Da:		Molecular weight to fit:	
populates these	230.10553	0.00026	230.10527	
boxes as you draw	Name:			
the structure.	N,N-Dimethyl-2-(2-methyl-1H-indol-3-yl)-2-oxoacetamide			
	ChemSpider	Apply FISh sco	oring Save Cancel	
	Opens	s the ChemSpider S	Search dialog box.	

Table 72 describes the parameters in the Compound Annotation Editor dialog box.

Parameter	Description
Description page	
Formula	Displays the elemental formula of the structure in the drawing area or the assigned formula.
Formula to Fit	Displays the elemental formula of the component found by the Find Expected Compounds node.
Molecular Weight/Error in Da	Displays the molecular weight (MW) of the structure in the drawing area and the difference between the structure's calculated MW and the MW for the selected compound—that is, the MW in the Molecular Weight to Fit box.
Molecular Weight to Fit	Displays the molecular weight (based on the formula) of the compound selected in the Compounds table or the Expected Compounds table.
Name	Displays the name of the compound from an online or local database search.
	To change the name, type alphanumeric text in this box.
Buttons and check box	at the bottom of the dialog box
ChemSpider	Opens ChemSpider Search dialog box for searching the ChemSpider database.
Apply FISh Scoring	Select this check box and click Save to run the FISh Scoring algorithm.
2	Saves the changes.
Save	

Table 72. Compound Annotation Editor parameters

Adding and Deleting Proposed Structures for a Compound

In a result file, every compound in the Compounds table or the Expected Compound table has a related Structure Proposals table.

To add or delete proposed structures for a compound, follow these topics:

- Adding Structure Proposals
- Deleting Structure Proposals

Adding Structure Proposals

- * To add structure proposals to a Structure Proposals table
- 1. Select a compound in the main Compounds table or the main Expected Compounds table.
- 2. Click Show Related Tables below the Compounds table or Expected Compounds table.
- 3. Click the Structure Proposals tab.
- 4. For each structure proposal, do the following:
 - a. Right-click anywhere below the Structure Proposals tab and choose **Structure Proposals > Add Structure Proposal**.
 - b. To edit the structure proposal, follow the instructions in "Editing Compound Annotations" on page 259.

Deleting Structure Proposals

* To delete a structure proposal from a Structure Proposals table

Right-click the row and choose **Structure Proposals > Delete Structure Proposal**.

Replacing an Annotation with a Structure Proposal

Annotations include the compound name, formula, annotation source, FISh coverage score, and structure.

To replace the annotations for a compound in the main result table

Right-click a row in the Structured Proposals table and choose **Structure Proposals > Use As Compound Annotation**.

Applying FISh Scoring

You can apply FISh scoring to a selected entry in the Compounds table, the Expected Compounds table, or the entries in the following compound-related search result tables: Structure Proposals, mzCloud Results, mzVault Results, Mass List Search Results, BioCyc Results, and ChemSpider Results. You can also apply FISh scoring from the Compound Annotation Editor dialog box.

The FISh scoring algorithm uses the structure in the structure column of the result table or the drawing area of the Compound Annotation Editor dialog box.

Follow these topics as needed:

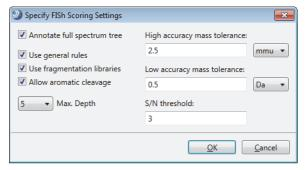
- Applying FISh Scoring from a Result Table
- Applying FISh Scoring from the Compound Annotation Editor Dialog Box
- Specifying the FISh Scoring Parameter Settings

Applying FISh Scoring from a Result Table

To apply FISh Scoring from a result table

- 1. Make sure that the entries of interest include structures.
- 2. To open the Specify FISh Scoring Settings dialog box (Figure 74), do one of the following:
 - To submit a compound in the Compounds or Expected Compounds table, right-click the selection and choose **Apply FISh Scoring**.
 - To submit a single selection in a Structure Proposals table, right-click the selection and choose **Structure Proposals > Apply FISh Scoring to Selection**.
 - To submit all of the entries in a Structure Proposals table, right-click the table and choose **Structure Proposals > Apply FISh Scoring to All**.
 - To submit an entry in a related search result table, right-click the table and choose Add to Structure Proposals and Apply FISh Scoring.

Figure 74. Specify FISh Scoring Settings dialog box



- 3. Set up the parameters (see "Specifying the FISh Scoring Parameter Settings" on page 264).
- 4. Click OK.

In the FISh Scoring Queue view to the left of the table, one job appears for each selected entry. For each entry that includes a structure, the application runs the FISh scoring algorithm. The run time increases as the complexity of the structure increases. When an entry does not include a structure, the job ends in failure and is highlighted with a red border.

Applying FISh Scoring from the Compound Annotation Editor Dialog Box

- * To apply FISh Scoring from the Compound Annotation Editor dialog box
- To open the FISh Scoring page of the Compound Annotations dialog box, right-click an entry in the Compounds table or the Expected Compounds table and choose Edit Compound Annotation.
- 2. Make sure that the drawing area on the Description page includes a structure.
- 3. Click the **FISh Scoring** tab.
- 4. Select the Apply FISh Scoring check box (Figure 75).

Figure 75. FISh Scoring page of the Compound Annotation Editor dialog box

Description FISh Scoring		
Annotate full spectrum tree	High accuracy mass tolerance:	
Use general rules	2.5	mmu 🔻
Use fragmentation libraries	Low accuracy mass tolerance:	
Allow aromatic cleavage	0.5	Da 🔻
5 🔹 Max. Depth	S/N threshold:	
	3	
ChemSpider	Apply FISh scoring Save	Cancel

- 5. Specify the parameter settings (see "Specifying the FISh Scoring Parameter Settings" on page 264).
- 6. Click Save.

Specifying the FISh Scoring Parameter Settings

- To set up the FISh scoring parameters
- 1. In the Specify FISh Scoring Settings dialog box or on the FISh Scoring page of the Compound Annotation Editor dialog box, make the following selections:
 - To annotate the full spectrum tree, select the **Annotate Full Spectrum Tree** check box.
 - To use the general fragmentation rules, select the Use General Rules check box.
 - To use the fragmentation libraries, select the Use Fragmentation Libraries check box.

Tip If time allows, select the **Use Fragmentation Libraries** check box. Using the fragmentation libraries provides significantly more structural information; however, it can also add a significant amount of processing time.

- To allow aromatic cleavage as one of the fragmentation steps, select the **Allow Aromatic Cleavage** check box.
- In the Max. Depth list, select the maximum number of steps allowed in the fragmentation pathway.
- 2. Use the default values or type new values in the following boxes:
 - For the FTMS scans, type a value in the High Accuracy Mass Tolerance box and select the appropriate units.
 - For the ITMS scans, type a value in the Low Accuracy Mass Tolerance box and select the appropriate units.
 - In the S/N Threshold box, type a value for the FTMS scans.

Table 73 describes the parameter settings for the FISh scoring algorithm.

Table 73. FISh scoring parameters

Parameter	Description		
Annotate Full Spectrum Tree	Annotates the full spectrum tree (MS/MS, MS ³ , and so on) in the Mass Spectrum view.		
Use General Rules	Uses the general fragmentation rules.		
Use Fragmentation Libraries	Uses the fragmentation libraries.		
Allow Aromatic Cleavage	Allows aromatic cleavage as one of the reaction steps.		
Max. Depth	Specifies the maximum number of fragmentation reactions to consider in the fragmentation pathway.		
	Default: 5; range: 1–20		
High Accuracy mass tolerance and units	Specifies the mass tolerance for FTMS data. Default: 2.5 mmu		
Low Accuracy Mass Tolerance and units	Specifies the mass tolerance for ITMS data. Default: 0.5 Da		
S/N Threshold	Specifies the signal-to-noise threshold for FTMS data. The FT mass analyzer calculates the S/N level for each centroid.		

Using Result Filters for Data Reduction and Creating Filter Sets

To show only the most pertinent data, use the Result Filters view to apply filters to the processed data. By default, the left pane of the Result Filters view lists the main tables in the current result file. The right pane displays the filters for the table that you select in the left pane.

Figure 76. Result Filters view with an empty filter for the Compounds table

② Result Filters	
ON Compounds ON Compounds per File ON Features ON MCOud Results ON Thiput Files	Compounds Add aroup Add property
Show all tables	Load Save Save As Clear All Clear Apply Filters

 Adds the visible related tables to the result filters list.

Note The filters for each table are independent of other table filters. For example, a retention time filter for the Merged Features table does not affect the entries in the other tables that include a retention time column. In addition, filtering only removes rows from the display; it does not update or change any of the calculated values.

To set up and apply the data reduction filters, see these topics:

- Setting Up, Applying, and Saving Filter Sets
- Creating a Result Filter for a Status Column
- Creating a Result Filter with an AND Logical Conjunction
- Creating a Result Filter with an OR Logical Conjunction
- Creating a Result Filter with Both of the Logical Conjunctions
- Loading a Saved Filter Set
- Result Filters View Parameters

Setting Up, Applying, and Saving Filter Sets

This topic describes how to set up, apply, and save a set of result filters (FILTERSET) for the result tables in a result file.

To set up, apply, and save a filter set

1. Open a result file.

2. From the application menu bar, choose View > Result Filters.

The Result Filters view opens as a floating window or as a docked view and displays the filter tree for the current table.

Each table has its own set of filter conditions.

3. (Optional) To display a filter tree that includes all the visible result tables in the result file, select the **Show All Tables** check box.

Note To change which tables are visible, use the Select Visible Tables dialog box.

4. In the Result Filters view, select the table of interest in the left pane.

The table name appears in the right pane. The following selection tree appears below the table name.

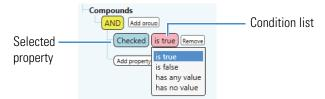
Table Name	
AND Add group	
Add property	Click to open a dropdown list of table columns.

5. Click Add Property.

A list that begins with the AND and OR logic selections followed by the columns in the selected table appears.

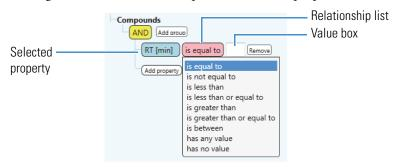
- 6. Select the table column (property) that you want to filter by.
- 7. Do any of the following:
 - For the Checked property, select a condition.

This figure shows the condition list for the Checked property.



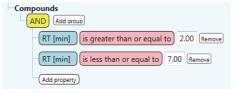
• For a numeric property, such as retention time (RT), select a mathematical relationship and type a value in the adjacent box, if applicable.

Tip When you select the Is Equal To relationship, type a numeric value to a minimum precision of two decimal places or a minimum precision that is equal to the number of decimal places that are displayed in the column, whichever is greater. For example, for any of the Area columns, type a numeric value with two decimal places, even though the Area column displays a numeric value with no decimal places.

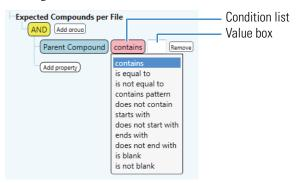


This figure shows the relationship list for numeric properties.

This figure shows a set of filter conditions that limits the displayed results to chromatographic peaks from 2 to 7 minutes.



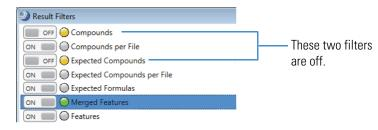
• For a non-numeric property, such as Parent Compound, select a condition and type a value in the value box if applicable.



This figure shows the condition list for non-numeric properties.

- 8. Do any of the following:
 - To apply the filter conditions in the current filter set, click **Apply Filters**.
 - To save a filter to a FILTERSET file, click **Save** or **Save As**. Then, browse to the location where you want to store the file, name the file, and click **Save**.
 - To turn off the filter conditions for a specific table, in the left pane of the Result Filters view, click **ON** to the left of the table name.

The indicator icon to the left of the table name turns from green to yellow, and the button displays OFF.



- To clear the filter conditions for a specific table, in the left pane of the Result Filters view, select the table. Then, click **Clear**.
- To clear all of the filters conditions in a filter set, click **Clear All**.

Creating a Result Filter for a Status Column

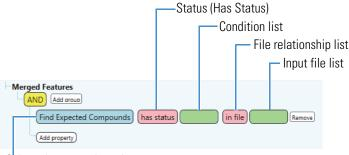
This topic describes how to set up a filter for a status column of a result table in a result file.

To set up filters for the status columns

- 1. In the Result Filters view, select one of these tables: Merged Features or Compounds.
- 2. Select a status property; for example, for the Merged Features table, select **Find Expected Compounds**.

From left to right, the following items appear:

- A (pink) list with one selection: Has Status
- A (green) condition list with no selection
- A (pink) file relationship list with the selection of In File
- A (green) input file list with no selection



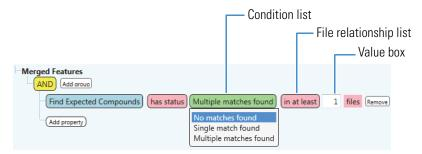
Selected status column in the Merged Features table

The input file list changes to a value box when you select anything other than In File in the file relationship list:

• If you select In File, select the file of interest in the input file list.

	Condition list	File relationship list
		Input file list
Merged Features		
Ion Conflict Status has status	Found only by 'Detect Unknown Compounds' node	in file Urine_0-4hr_01.raw (F3) Remove
(Add property)	No data to compare Not found by 'Detect Unknown Compounds' node Found only by 'Detect Unknown Compounds' node No conflict Conflicting ions Multiple ions per node	

• If you select a file relationship that requires additional input, a value box appears.



3. Select a status condition for the selected property (table column).

Ion Conflict Status	Find Expected Compounds, Detect Compounds, or Custom Explanations
No Data to Compare	No Matches Found
No Conflict	Single Match Found
Conflicting Ions	Multiple Matches Found
Multiple Ions Per Node	_
Not Found by Detect Unknown	_
Compounds node	
Found Only By Detect Unknown	_
Compounds node	

4. Select a file relationship and enter a value or select a file.

Creating a Result Filter with an AND Logical Conjunction

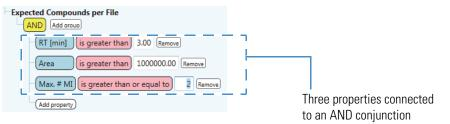
This topic describes how to use the AND logical conjunction in a result filter.

* To create a filter set using the AND logical conjunction

Note When you use the AND logical conjunction, all of the connected property conditions must be True.

- 1. In the Result Filter view, keep the **AND** logical conjunction as the first item in the filter tree.
- 2. For each property that you want to conjoin with the AND conjunction, click **Add Property**, select a property from the list, and set the property boundaries.

This figure shows a filter set that uses three properties conjoined with an AND conjunction. When you apply this filter set to the data in the Expected Compounds per File table, only those rows that meet all three conditions remain; that is, you see only those detected chromatographic peaks with a retention time greater than 3.00 minutes, with an integrated peak area greater than 1 000 000, and where at least two or more matching isotopes were found in the MS scans that make up the peak's data points.



Creating a Result Filter with an OR Logical Conjunction

This topic describes how to use the OR logical conjunction in a result filter.

* To create a filter set using the OR logical conjunction

Note When you use the OR logical conjunction, only one of the connected property conditions must be True.

- 1. Select the **OR** logical conjunction as the first item in the filter tree.
- 2. For each property that you want to conjoin with the OR conjunction, click **Add Property**, select a property from the list, and set up the property conditions.

This figure shows a filter set that uses two properties conjoined with an OR conjunction. When you apply this filter set to the data in the Expected Compounds table, those rows that meet at least one of the conditions remain—that is, you see the rows where the FISh coverage value is greater than or equal to 30 or where the FISh coverage column has no reported value.

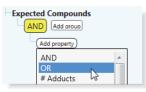


Creating a Result Filter with Both of the Logical Conjunctions

This topic describes how to use both of the logical conjunctions in a result filter.

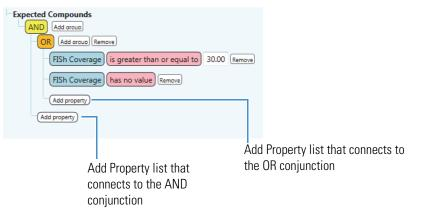
To create a filter set using both the AND and the OR logical conjunctions

- 1. Keep the AND logical conjunction as the first item in the filter tree.
- 2. To conjoin two properties with the OR conjunction, do the following:
 - a. Click Add Property and select OR from the dropdown list.



b. Set up the properties that you want to conjoin with the OR conjunction.

This example shows the condition where the FISh Coverage must be greater than or equal to 30 or have no value. (The FISh scoring algorithm does not calculate scores for DIA scans.)



3. For each property that you want to conjoin with the AND conjunction, click the **Add Property** list that connects to the AND conjunction, select a property from the list, and set up the property boundaries.

This figure shows a result filter that keeps chromatographic peaks that meet the following conditions:

- A FISh Coverage score that meets one of these conditions:
 - A FISh Coverage score that is greater than or equal to 30

-or-

No FISh Coverage score

-and-

• A Retention Time from 4 to 7 minutes

Expected Compounds
AND Add aroup
OR Add aroup Remove
FISh Coverage is greater than or equal to 30.00 Remove
FISh Coverage has no value Remove
Add property
RT [min] is greater than or equal to 4.00 Remove
RT [min] is less than or equal to 7.00 Remove
(Add property)

Loading a Saved Filter Set

This topic describes how to load the filters from a FILTERSET file.

To load a saved filter set

1. In the Result Filters view, click Load.

The Load Filter Set dialog box opens.

- 2. Browse to the appropriate folder and select the filter set of interest.
- 3. Click Open.

If the filter set contains filter conditions for tables that are not in the current result file, the application automatically hides the unused filters. You can modify the filter conditions for the applicable tables only.

Result Filters View Parameters

Use the Result Filters view to create data reduction filters for the tables in a result file. The Result Filters view is a floating window that can remain open while you work in other areas of the application.

Table 74 describes the panes, buttons, icons, and check box in the Result Filters view.

Table 74. Result Filters panes and buttons

Feature	Description
Left pane	Lists the main tables included in the current result file. An On/Off button and an indicator icon appear to the left of the table name.
ON/OFF button	Use to turn on or turn off the conditions for the associated main table.
Indicator icons	(⁽) Gray—Indicates that the table is unfiltered.
	(●) Green—Indicates that a filter has been applied to the table.
	() Yellow—Indicates that the table filter is off.

Right pane

Displays the filter settings for the selected table. You can modify these settings as described in "To set up, apply, and save a filter set" on page 266.

AND or OR	Specifies the logical connection between properties or groups.
Add Group	Adds a group.
Add Property	Adds a property.
Check box	
Show All Tables	Selecting this check box adds the related tables to the table list.
Buttons	
Load	Opens the Load Filter Set dialog box where you can select a saved filter set and open it.
Save	If a saved filter set is open, clicking Save overwrites the settings in the file with the current filter conditions in the Result Filters window.
Save As	Opens the Save Filter dialog box where you can define the file name and select a folder for a FILTERSET file.
Clear All	Clears all of the filter conditions for the current filter set.
Clear	Clears the current filter condition.
Apply Filters	Applies all of the filter conditions for the current filter set.

Viewing the Result Summary

In the Summaries view, you can view the following summaries:

- Workflow Summary
- Processing Messages Summary
- Filter Summary
- Study Summary
- Grouping & Ratios Summary

✤ To open the Summaries view

- 1. Open the result file of interest (see "To open a result file" on page 248).
- 2. From the menu bar, choose **View > Result Summary**.

The Summaries view includes these five pages: Workflow, Processing Messages, Filter, Study, and Grouping & Ratios (Figure 77).

Figure 77. Summaries view

② Summaries	
Workflow Processing Messages Filter Study Grouping & Ratios	
Search name: Target Omeprazole Search description: - Search date: 3/31/2017 4:53:48 PM Created with Discoverer version: 2.1	-
[Input Files (28)] >Select Spectra (29) >Create Analog Trace (36)	
[Select Spectra (29)] >Align Retention Times (30)	
[Align Retention Times (30)] >Find Expected Compounds (49)	
[Generate Expected Compounds (48)] >Find Expected Compounds (49)	
[Find Expected Compounds (49)] >Group Expected Compounds (50) >Merge Features (40)	
[Group Expected Compounds (50)] > Mark Background Compounds (45) > FISh Scoring (51)	
[Mark Background Compounds (45)]	
[FISh Scoring (51)]	
[Merge Features (40)]	
[Create Analog Trace (36)]	
[Differential Analysis (52)]	-
★	

Workflow Summary

To view the processing workflow used to create the active result file, open the Workflow page. This page lists the name of the processing workflow, the creation date for the result file (.cdResult), the raw data files (.raw) that were processed to create the result file, and the parameter settings for the workflow nodes.

For information about creating a processing workflow, see Chapter 6, "Creating and Editing Processing Workflows." For information about selecting one of the defined workflows, see "Nomenclature for the Defined Processing Workflows" on page 29.

Processing Messages Summary

To view a summary of the processes used to create the active result file, open the Processing Messages page.

Filter Summary

To view a summary of the filters used to reduce the data in the results window, open the Filter page. This page lists the name of the latest filter set (FILTERSET file type) that you applied to the result file and the filter conditions in the filter set. Use the Result Filters view to create filter sets (see "Using Result Filters for Data Reduction and Creating Filter Sets" on page 266).

Study Summary

To view a summary of the study settings for the input files that make up the result file, open the Study page. This page lists the following:

- Name and creation date of the study
- Directory location of the study
- Study factors and their values
- Sample names (Xcalibur RAW files) and their directory location

Grouping & Ratios Summary

To view a summary of the sample groups and ratios for the analysis, open the Grouping & Ratios page.

Figure 78 shows an example summary of the sample groups and ratios for an analysis. The summary lists the selected study variables, sample groups, and ratios in order from top to bottom.

Workflow Processing Messages Filter Study Grouping & Ratios Grouping:	*
Grouping: Study variable(s) for grouping: Time Points Sample Groups: n/a F1: Blank, n/a F2: Blank, n/a O-4hr	*
Study variable(s) for grouping: Time Points Sample Groups: n/a F1: Blank, n/a F2: Blank, n/a O-4hr	
Time Points Sample Groups: n/a F1: Blank, n/a F2: Blank, n/a 0-4hr	
Time Points Sample Groups: n/a F1: Blank, n/a F2: Blank, n/a 0-4hr	
Sample Groups: n/a F1: Blank, n/a F2: Blank, n/a 0-4hr	
Sample Groups: r/a F1: Blank, n/a F2: Blank, n/a 0-4hr	
n/a F1: Blank, n/a F2: Blank, n/a 0-4hr	
F1: Blank, n/a F2: Blank, n/a 0-4hr	
F1: Blank, n/a F2: Blank, n/a 0-4hr	
0-4hr	
F2: Sample 0.4hr	
F4: Sample, 0-4hr	
4-8hr	Ξ
F5: Sample, 4-8hr	
F6: Sample, 4-8hr	
8-12hr	
F7: Sample, 8-12hr	
F8: Sample, 8-12hr	
Predose	
F9: Sample, Predose F10: Sample, Predose	
rio: sample, riedose	
Ratios:	- 11
Katios:	
(0-4hr) / (Predose) (4-8hr) / (Predose)	
(8-12hr) / (Predose)	
4	

Figure 78. Grouping & Ratios Summary page

Exporting the Tabular Results to an External File

These topics describe how to export the contents of a result table to an external file:

- Shortcut Menu Commands for the Result Tables
- Exporting the Result Table Contents to a Spreadsheet
- Exporting the Result Table Contents to a Text File
- Exporting an Xcalibur Inclusion or Exclusion List
- Exporting the Contents of the Compounds Table to TraceFinder

Shortcut Menu Commands for the Result Tables

Table 75 describes the shortcut menu commands for the result tables.

Table 75.	Shortcut menu commar	ids for the res	sult tables ((Sheet 1 of 3)
-----------	----------------------	-----------------	---------------	----------------

Command	Description		
All main and related result tables			
Copy With Headers	Copies the current table row and its associated column headings to the Clipboard.		
Сору	Copies the current table row to the Clipboard. Does not copy the column headings.		
Clear Selection	Undoes any row selections. Clears the Chromatograms view, the Mass Spectrum view, or both of these views if they are populated with data.		
Cell Selection Mode	Turns on the cell selection mode. When the Cell Selection Mode is on, you cannot sort the table columns.		
Enable Column Fixing or Disable Column Fixing	Turns on the column pins. Pinning (or freezing) a column moves it to the left of the Checked column.		
Export > As Plain Text	Exports the data to a comma-separated values file.		
Export > As Excel	Exports the data to an Excel [™] spreadsheet file.		
All result tables with a Checked column			
Check Selected	Places a check in the selected row's check box.		
Check All	Selects the check boxes for all of the table rows.		
Uncheck Selected	Clears the selected row's check box.		
Uncheck All	Clears the check boxes for all of the table rows.		
Remove All Checkmarks in All Tables	Clears the check boxes in all the result tables.		
All result tables with expanding table headings			
Expand All Column Headers	Expands the collapsed column headings.		
Collapse All Column Headers	Collapses the expanded column headings.		

Command	Description	
Expected Compounds and Com	pounds tables	
Edit Compound Annotation	If the table does not already include the Name and FISh Coverage columns, adds these columns and opens the Compound Annotation Editor where you can name the compound, add a structure that matches the formula and molecular weight, run a ChemSpider search, and apply the FISh Scoring algorithm.	
Clear Compound Annotation	Removes the custom annotation.	
Export > As Xcalibur Inclusion/Exclusion List	Exports information about all of the compounds in the table or only the selected compounds in the table to a text file in the format required for the selected mass spectrometer. The information includes the m/z value of the monoisotopic ion and the start and stop times for the chromatographic peak.	
Export > Single Compound to an Existing mzVault Library	Exports a selected compound to an existing mzVault library.	
Export > Compounds to a New mzVault Library	Exports the selected compounds to a new mzVault library.	
Compounds table		
Export > Export to TraceFinder	Exports information about all of the compounds, only the named compounds, or only the checked compounds in the table to a CSV file in a format appropriate for the TraceFinder application.	
Export > As Mass List	Exports the selected items to a new mass list.	
Add Selected Compounds to Existing Mass List	Adds the selected compounds to the specified mass list.	
Molecular Networks (Beta)	Opens the "molecular networking" viewer in your default web browser.	
Structure Proposals table (rela	ted to the Compounds and Expected Compounds tables)	
Structure Proposals > Add Structure Proposal	Adds a new row to the Structure Proposals table that includes the formula and molecular weight from the selected row in the main Compounds or Expected Compounds table.	
	You can type a name in the Name column and a description in the Comments column. Double-click the new row to open the Compound's Annotation Editor where you can draw the compound's structure, open a structure file, or run a ChemSpider search.	

 Table 75.
 Shortcut menu commands for the result tables (Sheet 2 of 3)

Command	Description	
Structure Proposals > Edit Structure Proposal	Opens the Compound Annotation Editor where you can name the compound, add a structure that matches the formula and molecular weight, run a ChemSpider search, and apply the FISh Scoring algorithm.	
Structure Proposals > Delete Structure Proposal	Removes the selected row from the Structure Proposals table.	
Structure Proposals > Use as Compound Annotation	Replaces the annotations in the selected row of the Compounds or Expected Compounds table with the annotations in the current row of the Structure Proposals table.	
Structure Proposals > Apply FISh Scoring to Selected	Opens a dialog box where you can specify the settings for the FISh scoring algorithm and submit the information in the selected row.	
	To apply FISh scoring, the selected row in the Structure Proposals table must include a formula or a structure.	
Structure Proposals > Apply FISh Scoring to All	Opens a dialog box where you can specify the settings for the FISh scoring algorithm and submit all the rows in the Structure Proposals table.	
All related tables that have a c	orresponding main table	
Go to Same Item in Main Table	Opens the main table with the same name as the related table and selects the corresponding table row in the main table.	

Table 75. Shortcut menu commands for the result tables (Sheet 3 of 3)

Related Topics

- Copying Table Entries to the Clipboard
- Freezing Table Columns
- Exporting the Result Table Contents to a Text File
- Exporting the Result Table Contents to a Spreadsheet
- Exporting an Xcalibur Inclusion or Exclusion List

Exporting the Result Table Contents to a Spreadsheet

You can export the contents of any of the result tables to a spreadsheet file.

* To export the contents of a result table to a spreadsheet file

1. Right-click the result table that you want to export and choose **Export > Export to Excel**.

The Export to Excel dialog box opens (Figure 79). The default storage path is C:\Users\Public\Documents*result file name*. The default selection in the Level 1 box is the name of the active result table.

Figure 79. Export to Excel dialog box with the Compounds table selected

Export to Exce	l		—
Path C:\Users\Public	\Documents\Unknown Expected De	scriptive Statist	tics.xlsx
	iated tables to be exported	• •	Options Checked items only Open file after export
		Ex	port Cancel

2. If necessary, select a different result table from the Level 1 list.

The list includes all the main tables in the result file.

3. Click the browse icon next to the Path box.

A second Export to Excel dialog box (where you specify the location and file name for the file) opens in front of the first Export to Excel dialog box.

Export to Excel			
○○ - ○ < M	Documents CD_Exports	✓ ✓ Search CD_Exports	٩
File <u>n</u> ame:	Expected Compounds per File.xlsx		•
Save as <u>t</u> ype:	XLS (Excel Spreadsheet) (*.xlsx)		-
<u> B</u> rowse Folders ■ Compare Folders ■ Compare Folders ■ Compare Folders ■ Compare Folders Compare Folders		Save Cancel	

4. Browse to the folder where you want to store the file.

By default, the File Name box displays the name of the result table that you are exporting.

- 5. Overwrite or rename the file as appropriate, select the spreadsheet type (**XLS** or **Microsoft Excel File**), and click **Save**.
- 6. In the Export to Excel dialog box (for selecting the table rows to be exported), select related tables from the Level 2 and Level 3 lists as appropriate.

- 7. In the Options area, make the following selections:
 - To export only those rows that are checked in the result table or tables, select the **Checked Items Only** check box. Otherwise, leave the check box clear to export all of the table rows.
 - To automatically open the newly created spreadsheet after you export it, select the **Open File After Export** check box.
- 8. Click **Export** to export the data to a spreadsheet.

When the export is finished, you receive a confirmation message with the file's location.

9. At the prompt, click **OK**.

Exporting the Result Table Contents to a Text File

You can export the contents of any of the result tables to a text file.

* To export the contents of a result table to a text file

 Right-click the result table that you want to export and choose Export > Export to Text File.

The Export to CSV File dialog box opens. The File Name box displays the name of the selected result table.

File <u>n</u> ame:	Compounds.csv	,
Save as <u>t</u> ype:	CSV (Comma delimited) (*.csv)	,

2. Select the folder where you want to store the file, name the file as necessary, and click **Save**.

The text file (.csv) appears in the selected folder.

Exporting an Xcalibur Inclusion or Exclusion List

You can export the contents of the Compounds table or the Expected Compounds table to an Xcalibur Inclusion/Exclusion list, and then import this list into an instrument method that controls your Thermo Scientific mass spectrometer.

To export an Xcalibur inclusion or exclusion list

Right-click the Compounds table or the Expected Compounds table and choose Export
 > As Xcalibur Inclusion/Exclusion List.

The Export Xcalibur Inclusion/Exclusion List dialog box opens (Figure 80).

C·\LIsers\Public\Documents\OC	Normalization 50_30_2_million.txt
Options Left RT tolerance [min]: 1 Right RT tolerance [min]: 1 Checked items only Include isotopic peaks	Instrument Instru
LTQ Orbitrap Options Mass precision (decimals): 5 Max. concurrent entries: 5 Remove charge	10

Figure 80. Export Xcalibur Inclusion/Exclusion List dialog box

- 2. Make the appropriate selections and entries (Table 76).
- 3. Click Export.

The application attempts to save the text file to the specified location. If the file name conflicts with an existing file, a confirmation message appears. In the absence of a conflicting file, a completion message appears indicating that the exported file is in the selected folder.

- 4. If a confirmation message appears, do one of the following:
 - To overwrite the existing file, click **Yes**, and then click **OK** when the completion message appears.
 - To cancel the export, click No.

Table 76 describes the parameters in the Export to Xcalibur Exclusion List dialog box.

	Table 70. Export to Adalbar inclusion/Exclusion List dialog box parameters (Sheet 1 of S)		
Parameter	Description		
Path	Specifies the file name and directory path of the text file that contains the inclusion/exclusion list for your Xcalibur instrument method.		

Table 76. Export to Xcalibur Inclusion/Exclusion List dialog box parameters (Sheet 1 of 3)

Table 76.	Export to Xcalibur	Inclusion/Exclusion List of	dialog box parameter	s (Sheet 2 of 3)

Parameter	Description		
Options			
IMPORTANT An Orbitrap Fusion mass spectrometer accepts a retention time range of 0.01 to 999 minutes.			
	Make sure that the retention time window for each compound falls within the retention time range of the instrument method.		
• Expected RT – Let method	ft RT Tolerance > Minimum retention time for the instrument		
• Expected RT + Rig method	ght RT Tolerance < Maximum retention time for the instrument		
Left RT Tolerance [min]	Specifies the minimum start time for the chromatographic peak. If the Expected RT minus the Left RT Tolerance setting is less than zero, the application exports a value of zero.		
	Default: 1 Range: 0.001 to 1000		
Right RT Tolerance	Specifies the maximum stop time the chromatographic peak.		
[min]	There is no error checking for the calculated maximum retention time.		
	Default: 1 Range: 0.001 to 1000		
Checked Items Only	Specifies that the application exports only the selected compounds to the named text file.		
	Default: Clear		
Include Isotopic Peaks	Adds an entry for each isotopic spectral peak.		
LTQ Orbitrap Options			
Mass Precision	Specifies the mass precision.		
(Decimals)	For the LTQ Orbitrap, the mass precision of the exported data must match the required mass precision for your Xcalibur instrument methods. You specify the required mass precision for Xcalibur instrument methods in the Instrument Configuration dialog box of the Foundation platform.		
	Default: 5; range: 0 to 5		

Parameter	Description	
Max. Concurrent Entries	Specifies the maximum number of entries with overlapping the windows.	
	Default: 500; range: 1 to 2000	
Remove ChargeSpecifies whether the application exports the <i>m/z</i> valion for each detected compound or the neutral mass detected compound.		
	Default: Clear	
Instrument		
LTQ Orbitrap	Exports the list in the appropriate format for the LTQ Orbitrap instrument control software.	
Q Exactive	Exports the list in the appropriate format for the Q Exactive instrument control software.	
	The mass list includes the formula of each compound in the Comment column when the formula is available.	
Orbitrap Fusion	Exports the list in the appropriate format for the Orbitrap Fusion instrument control software.	
	If the table contains both positive and negative scans, the application creates two mass lists.	
	The maximum number of target compounds for an Orbitrap Fusion mass list is 50 000. If the table includes more than 50 000 compounds, filter the table or check the compounds of interest before you export the mass list.	
Buttons and check box		
Open File(s) after Export	Specifies that the file opens after the application completes the export.	
Export	Exports the specified information to a text file.	
Close	Closes the dialog box.	

Table 76. Export to Xcalibur Inclusion/Exclusion List dialog box parameters (Sheet 3 of 3)

Exporting the Contents of the Compounds Table to TraceFinder

You can export the contents of the Compounds table to a CSV file that the TraceFinder application can use for data processing.

* To export a compounds list for the TraceFinder application

1. Right-click the Compounds table and choose Export > As TraceFinder List.

The Export to TraceFinder dialog box opens. The default folder is either *drive*:\Users\Public\Documents or the last folder you selected. The default file name is the name of the active result file.



② Export to TraceFinder	X
Path:	
C:\Users\Public\Documents\Pattern	Scoring and compound class 🛄
Options	
Exclude items without name	
Checked items only	
Open file after export	
	Export Cancel

- 2. Select the folder where you want to store the file.
- 3. Do the following as needed:
 - To exclude unnamed compounds, select the **Exclude Items Without Name** check box.
 - To include only the checked compounds, select the Checked Items Only check box.
 - To automatically display the exported compounds list, select the **Open File After Export** check box.
- 4. Click Export.

A status box appears when the export process finishes.

5. Click OK.

Exporting Spectral Data to a New or Existing mzVault Library

To export spectral data to a new or existing mzVault Library file, see these topics:

- Adding a Compound to an Existing mzVault Library
- Creating a New mzVault Library

Adding a Compound to an Existing mzVault Library

From a Compounds or Expected Compounds result table, use the Export > Single Compound to an Existing mzVault Library shortcut menu command to export a selected compound to an existing mzVault library.

Tip The application comes with an empty library named Custom mzVault Library.db. You can create other libraries by clicking New on the Libraries > Spectral Libraries page.

- To add a compound to an existing mzVault Library
- 1. Open a result file (see "Opening, Closing, and Updating Result Files" on page 248).
- 2. Click the **Compounds** tab or the **Expected Compounds** tab.
- 3. Right-click a compound in the table and choose **Export > Single Compound to an Existing mzVault Library**.

The Export to mzVault Library dialog box opens (Figure 82). Its Spectra view displays the available fragmentation spectra for the selected compound.

) Export to mzVa	ult Library									×
Compound L-(-)-Methionine C5 H11 N O2 S		Spectra	a (!:, #967, RT=1.52	8 min, MS2,	FTMS (🔊	700	_ddMS2 (F2) #967, RT=1.52	8 min, MS2, F1 172.04037	FM: •
_S	OH VH ₂	*	111		Þ	 Intensity [counts] (10,3) 100 100<!--</th--><th>66.03 109:05 60 80</th><th>011 172.09</th><th></th><th>→ 4 III</th>	66.03 109:05 60 80	011 172.09		→ 4 III
mzVault Library Selected library: Custom mzVault Library.db										
Structure	Name	Formula	Molecular Weight	Best Match	mzVault II	D mzClo	oud ID	ChemSpider ID	KEGG ID	
<u>A</u> a •	<u>A</u> a •	<u>A</u> a ▼		= -	= -	=	•	= -	<u>A</u> a •	
						Add as Nev	v (Add to Selected	Clos	ie 🛛

Figure 82. Export to mzVault Library dialog box populated with a Compounds table compound

- 4. Do the following:
 - In the spectrum tree, select the check boxes for the spectra that you want to add to the compound entry.
 - In the mzVault Library area, select the existing mzVault library from the Selected Library list.

- 5. Do one of the following:
 - To add the spectra to an existing compound entry in the library, select the entry in the Similar Compounds Found in Library list. Then, click **Add to Selected**.
 - To add the compound and the selected spectra as a new compound entry, click **Add as New**.
- 6. Close the dialog box.

Creating a New mzVault Library

From a Compounds or Expected Compounds result table, use the Export > Compounds to a New mzVault Library shortcut menu command to export all or selected compounds to new mzVault library.

- * To export compounds to a new mzVault library
- 1. Open a result file (see "Opening, Closing, and Updating Result Files" on page 248).
- 2. Click the **Compounds** tab or the **Expected Compounds** tab.
- 3. (Optional) Filter the table to display only the compounds of interest or select the check boxes for the compounds of interest.
- 4. Right-click the table and choose **Export > Compounds to a New mzVault Library**.

The New mzVault Library dialog box opens. By default, the application populates the Library Name box with the name of the opened result file.

Export to New mzVault Library	
Library Name:	
Target Omeprazole.db	
Options	
Exclude items without name	
Checked items only	
	Export Cancel

- 5. Do the following:
 - Type a new library name or keep the default name.
 - In the Options area, select whether to exclude compounds without a name, export only the compounds with selected check boxes, or both.
- 6. Click Export.
- 7. At the prompt, click **OK**.

The new library appears in the Spectral Libraries list.

8. To modify a library with the mzVault 2.2 application, export it to another folder. Then, use the Replace command to replace the existing library with the modified library.

Exporting Compounds to a New or Existing Mass List

You can create new mass lists or add compounds to existing mass lists by exporting compounds from the Compounds table of a result file.

Follow these topics as needed:

- Exporting Compounds from the Compounds Table to a New Mass List
- Exporting Compounds from the Compounds Table to an Existing Mass List

Exporting Compounds from the Compounds Table to a New Mass List

- To create a new mass list by exporting a set of compounds from the Compounds table
- 1. Open the result file of interest.
- 2. Open the Compounds table.
- 3. (Optional) To export only the checked compounds, select the check boxes for the compounds of interest in the Checked column.
- 4. Right-click the Compounds table and choose **Export > As Mass List**.
- 5. In the Export to New Mass List dialog box, do the following:
 - a. In the Mass List Name box, type a name for the mass list.
 - b. In the Options area, do any of the following:
 - To export only checked compounds, select the Checked Items Only check box.
 - To export only named compounds, select the **Exclude Items without Name** check box.
 - To include the retention time information in the mass list, select the **Export Retention Time** check box.
 - c. Click Export.
- 6. At the prompt, click **OK**.

The new mass list appears in your Mass Lists library.

Exporting Compounds from the Compounds Table to an Existing Mass List

To export compounds to an existing mass list

- 1. Open a result file of interest.
- 2. Open a main or related Compounds table.
- 3. Highlight the compounds of interest by using the SHIFT and CTRL keys.

4. Right-click your selection and choose Export > Add Selected Compounds to Mass List.

The Export to Existing Mass List dialog box opens.

Figure 83. Export to Existing Mass List dialog box

Export to Existing Mass List						
Mass list name:						
	•					
Options						
 Do not add duplicate names Export retention time 						
Entries Matching						
Mass tolerance: 1.0	ppm					
RT tolerance: 0.1	min					
(Export Cancel					

- 5. Do the following:
 - In the dropdown Mass List Name list, select the mass list where you want to add the selected compounds.
 - (Optional) To add the retention time information to the mass list, select the **Export Retention Time** check box.
- 6. To avoid exporting named compounds that are already in the mass list, do one of the following:
 - To exclude any named compound with a name that matches that of a named compound in the existing mass list, select the **Do Not Add Duplicate Names** check box.

Note This feature is case sensitive.

-or-

a. To exclude a named compound only if its mass and retention time match the duplicate compound in the mass list within specified tolerances, clear the **Do Not Add Duplicate Names** check box.

The Mass Tolerance and RT Tolerance boxes become available.

b. Enter the tolerances that you want the application to use to exclude duplicate named compounds from being exported to the mass list.

Note If you do not want to exclude a named compound unless both its mass and retention time are an exact match to the named compound in the mass list, set both tolerance values to 0.

Copying or Saving Graphical Views for Publication

You can copy the contents of a graphical view (see "Working with the Graphical Result Views" on page 293) to the Clipboard as editable text or as an image, or to an external file.

- To save a graphical view as a raster image, save it as one of these file types: PNG, GIF, JPG, TIF, or BMP.
- To save a graphical view as a vector image, save it as an EMF file.

Table 77 lists the shortcut menu commands for copying data to the Clipboard or an external file.

Table 77.	Commands for	copying an ima	age of a graphical	view (Sheet 1 of 2)

Graphical view	Copy the image to the Clipboard	Copy the data points to the Clipboard	Copy the data to an image file	Copy the data to a TEXT or CSV file
Chromatograms view	Copy > Image	Copy > Points	Export > Image As	Export > Points As
Mass Spectrum view	Copy > Image	Copy > Points—Copies the scan label and the <i>m/z</i> and intensity values for annotated centroids to the Clipboard.	Export > Image As	Export > Points As
		Copy > Raw Points—Copies the m/z and intensity values for all centroids to the Clipboard. Does not copy the scan label.		
Scatter Chart	Copy > Image	Copy > Data	Export > Image As	Export > Data As
Histogram Chart	Copy > Image	Copy > Data	Export > Image As	Export > Data As
Bar Chart	Copy > Image	Copy > Data	Export > Image As	Export > Data As
Pie Chart	Copy > Image	Copy > Data	Export > Image As	Export > Data As
Trend Chart	Copy > Image	Copy Information to Clipboard	Export > Image As	Export > Data As
Principal Component Analysis view	Copy > Image	Copy > Points	Export > Image As	Export > Points As (CSV file)
PLS-DA view	Copy > Image	Copy > Points	Export > Image As	Export > Points As
		Copy > Point Details		(CSV file)
		**		Export > Point Details As
Descriptive Statistics view	Copy > Image	Copy Information to Clipboard	Export > Image As	Save Information As (CSV file)

Graphical view	Copy the image to the Clipboard	Copy the data points to the Clipboard	Copy the data to an image file	Copy the data to a TEXT or CSV file
Differential Analysis view	Copy > Image	Copy > Points Copy > Point Details (Use this command if the corresponding table includes structures.)	Export > Image As	Export > Points As Export > Point Details As
KEGG Pathways	Сору	N/A	Save Picture As (a PNG or BMP file)	N/A
Retention Time Correction	Copy > Image	Copy > Points	Export > Image As	Export > Points As
Compound Area Corrections	Copy > Image	Copy > Points	Export > Image As	Export > Points As

Table 77. Commands for copying an image of a graphical view (Sheet 2 of 2)

Working with the Graphical Result Views

To work with the graphical views that are available when a result page is active, see these topics.

Contents

- Working with the Chromatograms View
- Working with the Mass Spectrum View
- Working with the Result Charts
- Plotting the Data as a Histogram Chart
- Working with the Trend Chart View
- Working with the Mass Defect Plot View
- Working with the Principal Component Analysis View
- Working with the Descriptive Statistics View
- Working with the Differential Analysis View
- Working with the Partial Least Squares Discriminant Analysis View
- Viewing KEGG Pathways
- Viewing BioCyc Pathways
- Viewing Metabolika Pathways
- Viewing the Corrected Retention Times of the Alignment Features
- Viewing the QC-Based Compound Area Correction Curves
- Viewing the FISh Scoring Queue
- Running a Hierarchical Clustering Analysis
- Running an mzLogic Analysis
- Working with the Isotopologues Distribution Chart

For information about using the Result Filters view, see "Using Result Filters for Data Reduction and Creating Filter Sets" on page 266. For information about viewing a summary of the analysis parameters for a result file, see "Viewing the Result Summary" on page 275.

Working with the Chromatograms View

When you initially open a result file, the Chromatograms view displays the XIC traces for the compound in the first row of the Compounds table or the Expected Compounds table. By default, the display zooms in on the detected peaks for the selected rows.

The Chromatograms view consists of a collapsible pane on the left and the graphical view on the right. Right-clicking the graphical view opens a shortcut menu (see Table 79).

Follow these topics as needed:

- Viewing a Chromatogram
- Adding a Chromatogram Plot
- Overlaying Chromatogram Plots
- Changing the Grouping in the Collapsible Pane
- Hiding the Traces for a Study Variable Value
- Updating All the Chromatogram Plots Simultaneously
- Manually Integrating Chromatographic Peaks
- Chromatograms View Shortcut Menu Commands

Viewing a Chromatogram

You can display chromatogram traces (a plot of intensity versus time) by selecting a row in any of these tables: Compounds, Compounds per File, Expected Compounds, Expected Compounds per File, Expected Formulas, Merged Features, Expected Features, FISh Trace Fragments, Specialized Traces, or Manual Peaks.

To view a chromatogram

- 1. Open a result file that contains chromatographic data.
- 2. If the Chromatograms view is closed, open it by choosing **View > Chromatogram** from the menu bar.

Note When you select a row for a filtered study file, the Chromatograms view does not display a trace. For example, if the Blank check box is clear under Filter By Sample Type and you select a row for a Blank sample, the Chromatograms view remains empty. By default, the Blank Sample Type check box is clear until you select it.

- 3. Do one of the following:
 - Click the **Compounds** tab or the **Expected Compounds** tab and select a row.

The Chromatograms view displays overlaid traces of the chromatographic peaks that were detected across the input files for the molecular weight and retention time (MW \times RT dimensions) listed in the selected table row. Each trace is a composite of the adducts found. For the Expected Compounds table, the chromatographic peak for each row is also derived from the same parent compound and reaction steps.

• Click the **Expected Formulas** tab and select a row.

The Chromatograms view displays overlaid traces for all of the chromatographic peaks found for an expected elemental composition (same MW, parent compound, and elemental composition).

• Click the Merged Features tab and select a row.

The Chromatograms view displays overlaid traces of the chromatographic peaks detected by the Detect Compounds and Find Expected Compounds nodes for the selected feature (same $m/z \times RT$ dimensions).

• Click the Features tab or the Expected Features tab and select a row.

The Chromatograms view displays the integrated chromatographic peak for the selected table row ($m/z \times RT$ dimensions).

• Click the **Compounds per File** tab or the **Expected Compounds per File** tab and select a row.

The Chromatograms view displays the integrated chromatographic peak for the selected table row (MW \times RT dimensions). Each trace is a composite of its related adducts, and the peak area for a compound is the sum of the areas for its related adducts (parent area for the adducts listed in the Expected Features table or the Features table).

• Click the FISh Trace Fragments tab and select a row.

The Chromatograms view displays a trace for the structure displayed in the selected table row. The FISh Trace Fragments table appears in the main table set when you select True for Individual Traces in the Create FISh Trace node of a processing workflow.

• Click the Specialized Traces tab and select a row.

Note The Specialized Traces table contains traces generated by these nodes: Create Mass Trace, Create Analog Trace, Create Pattern Trace, and Create FISh Trace.

Workflow node	Generates any of these trace	
Create Mass Trace	An extracted ion (mass range) chromatogram (XIC), a base peak chromatogram (BPC), or a total ion chromatogram (TIC)	
Create Analog Trace	A UV-Vis trace from a UV-Vis or PDA detector, up to three traces from a PDA detector, or an analog trace from an LC detector that you connected to one of the analog input channels of a Thermo Scientific mass spectrometer	
Create Pattern Trace	A TIC trace of the summed intensities of the mass spectral peaks (across the entire scan) that match the user-defined isotope pattern	
Create FISh Trace	• A summed FISh trace of all the matching fragment ion scans [data-dependent acquisition (DDA) or da independent acquisition (DIA)] when you select Tru in the Summed Trace list.	
	• An individual FISh trace for each fragment ion when you select True in the Individual Traces list. To view the individual trace for each fragment ion, see the FISh Trace Fragments table.	

Table 78. Wo	rkflow nodes	that generate s	pecialized traces
--------------	--------------	-----------------	-------------------

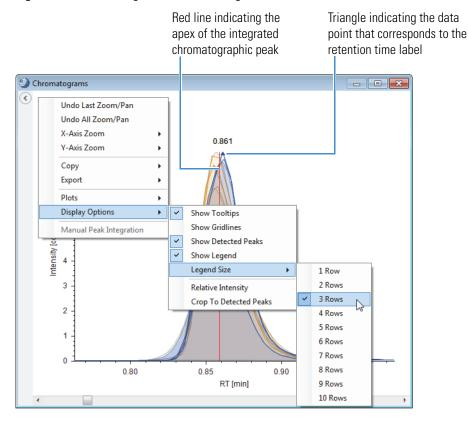
• Open the related tables, click the Filled Gaps tab, and select a row.

Tip By default, the Filled Gaps table is hidden. To display this table, open the Select Visible Tables dialog box, select the **Filled Gaps** check box, and click **OK**. For more information, see "Filled Gaps Table" on page 411.

- 4. To determine the origin of a trace in a result file that includes multiple input files, right-click the Chromatograms view and choose **Display Options > Show Legend**.
- To decrease or increase the number of legends displayed, right-click the Chromatograms view and choose Display Options > Legend Size > #Rows, where # is an integer value from 1 to 10 (Figure 84).

In Figure 84, the vertical red line indicates the peak apex of an integrated chromatographic peak. The triangle below the retention time label indicates the corresponding data point in the XIC trace.





When the number of legend lines becomes too large for the available display space, the application displays an empty view with the following text:

Not enough space for drawing the chart properly.

Adding a Chromatogram Plot

To add a plot to the Chromatograms view

- 1. Open a result file that contains chromatographic data.
- 2. Right-click the Chromatograms view and choose Plots > Add Plot.

The new plot appears below the original plot. Only the screen size of your computer monitor limits the maximum number of plots that you can add to the Chromatograms view. When you exceed the maximum number of plots, the plots disappear from the view.

In a Chromatograms view with more than one plot, a light blue bar on the left border highlights the active plot (Figure 85). If you right-click the Chromatograms view and choose Plots > Remove Plot, the application removes the active plot.

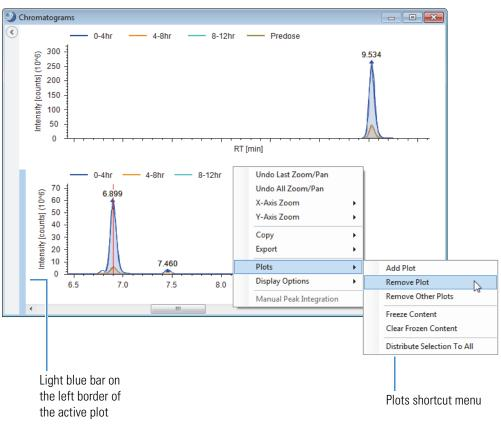


Figure 85. Chromatograms view with two plots

Overlaying Chromatogram Plots

* To overlay multiple chromatograms in one chromatogram plot

- 1. Open a result file that contains chromatographic data.
- 2. Do one of the following:
 - Hold down the SHIFT key and select a range of contiguous rows.
 - Hold down the CTRL key and select contiguous or noncontiguous rows one by one.
 - Hold down the SHIFT key and press the Down Arrow on the keyboard.

Tip Take care to avoid clicking an editable column in the result table, as doing so undoes the row selection and sets the focus to the table cell.

Changing the Grouping in the Collapsible Pane

To change the grouping of the chromatogram traces

- 1. Open the collapsible left pane by clicking the icon,
- 2. Under Group By, select or clear one or more of the check boxes.
- 3. To display all of the study file traces in different colors, clear all of the check boxes under Group By, or under Group By, select only the **Samples** check box.

Hiding the Traces for a Study Variable Value

***** To hide the chromatogram trace or traces for a study variable value

Under Filter By, clear the check box for the study variable value.

Updating All the Chromatogram Plots Simultaneously

- To update all the plots in the Chromatograms view simultaneously
- 1. Open a result file that contains chromatographic data.
- 2. Add two or more plots to the Chromatograms view.
- 3. Right-click the view and choose **Plots > Distribute to All Selections**.

As you select different table rows, all of the plots update. When Distribute to All Selections is not enabled, only the active plot updates.

Manually Integrating Chromatographic Peaks

- To manually integrate chromatographic peaks
- 1. Open a result file that contains specialized traces.

Note These workflow nodes generate specialized traces: Create Mass Trace, Create Analog Trace, Create Pattern Trace, and Create FISh Trace.

- 2. In the Specialized Traces table, select the trace of interest.
- 3. Right-click the Chromatograms view and choose Manual Peak Integration.

Two red dashed lines appear and the integrated peak area appears in blue.

4. Drag one line to the beginning of the chromatographic peak and the other line to the end of the chromatographic peak.

The integrated peak area changes as you move the start and end points of the chromatographic peak.

5. Place the cross-hair cursor on the peak.

A pop-up box appears with the selected retention time range and the Apply and Cancel buttons (Figure 86).

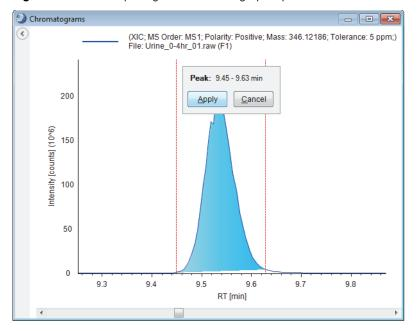


Figure 86. Manually integrated chromatographic peak

6. To add the manual peak to the Manual Peaks table, either click **Apply**, or on the keyboard, press A.

If the result file did not already contain a Manual Peaks table, the new table appears in the main table tab set. The Manual Peaks table contains a row for the new manual peak.

Chromatograms View Shortcut Menu Commands

Table 79 describes the shortcut menu commands for the Chromatograms view.

Table 79. Shortcut menu commands for the Chromatograms view (Sheet 1 of 4)

Command	Description	
Note The default zoom settings are as follows:		
• X-axis Zoom > To Detected Peaks (Active Plot)		
• Y-axis Zoom > Auto Scale Y-axis		
Undo Last Zoom/Pan	Undoes the last zoom or pan movement.	
Undo All Zoom/Pan	Zooms out to the full data acquisition time for the chromatogram on the x axis and the height of the largest chromatographic peak on the y axis.	

Description Maintains the same x-axis zoom range as you select different table rows. Overrides the Zoom to Detected Peaks command. Displays the full data acquisition time for the chromatogram. Zooms the x axis to the detected peaks for the selected table rows. Zooms the x axis to display the detected peaks in all of the plots.
rows. Overrides the Zoom to Detected Peaks command. Displays the full data acquisition time for the chromatogram. Zooms the <i>x</i> axis to the detected peaks for the selected table rows.
Zooms the <i>x</i> axis to the detected peaks for the selected table rows.
-
Zooms the x axis to display the detected peaks in all of the plots.
Default selection—Scales the y axis to the maximum intensity within the current <i>x</i> -axis (retention time) zoom range.
Manual zooming of the <i>y</i> axis is unavailable in this mode.
Scales the <i>y</i> axes of all the plots to the same scale. Manual zooming on both axes is available.
To reset the <i>y</i> -axis scaling, choose Y-axis Zoom > Auto Scale Y-Axis , and then choose Undo All Pan/Zoom .
Only changes the <i>y</i> -axis scaling of the current plot. Manual zooming on both axes is available.
Copies an image of the Chromatograms view (including the legend) to the Clipboard.
You can paste the image into a Microsoft Office document as a raster image or into a vector-drawing program as a vector image.
Copies the data as a two-column list of data points and copies the scan header. The first column lists the retention time and the second column lists either the relative intensity or the counts.
Opens the Save As dialog box where you can specify a file name and save the contents of the Chromatograms view as one of these selectable image formats: EMF, PNG, GIF, JPG, TIFF, or BMP. The EMF format is a vector image.

 Table 79.
 Shortcut menu commands for the Chromatograms view (Sheet 2 of 4)

Command	Description		
Exports > Points As	Opens the Save As dialog box where you can specify a file name and save the contents of the Chromatograms view as a text file. The default file name is Chart.txt.		
	When the chromatogram is a plot of relative intensity versus retention time, the application saves the data points as a two-column list. The first column lists the retention time and the second column lists the relative intensity (%). When the chromatogram is a plot of area (in counts) versus retention time, the second column lists the area for the peaks.		
Plots > Add Plot	Adds an empty, active plot to the Chromatograms view. Only the screen size limits the maximum number of displayed plots. When you reach the screen's limit, the Chromatograms view appears to be empty and the following message appears:		
	Not enough space for drawing chart properly.		
Plots > Remove Plot	Removes the active plot, which has a gray border.		
	Adding more than one plot activates this command.		
Plots > Remove Other Plots	Removes all of the plots in the Chromatograms view.		
	Adding more than one plot to the Chromatograms view activates this command.		
Plots > Freeze Content	Keeps the chromatogram of the currently selected row in the view when you select another row (in the current table or another result table). The application uses dashes to distinguish the frozen chromatogram trace.		
	Frozen peak with dashes		
	Chromatograms		
	F4: Urine_0-4hr_02 F3: Urine_0-4hr_01		
	Clears the frozen chromatogram from the view.		

Table 79. Shortcut menu commands for the Chromatograms view (Sheet 3 of 4)

Content

Command	Description	
Plots > Distribute Selection to All	Updates all of the plots simultaneously as you select different table rows. When Plots > Distribute Selection to All is not enabled, only the active plot updates.	
Display Options > Show Tooltips	 Displays a pop-up box with the following information, from top to bottom, when you place the cursor, +, over a chromatographic peak: Parent compound and any applicable transformations Chemical formula of the peak component Molecular weight of the peak component Selected retention time, in minutes Intensity, in counts, of the selected point on the chromatogram trace File name of the input file 	
Display Options > Show Gridlines	Adds grid lines to the Chromatograms view.	
Display Options > Show Detected Peaks	Uses a fill color for the integrated area under the detected chromatographic peaks.	
	Turning off this command removes the fill color.	
	Tip To see the peaks underneath the larger peaks in a set of overlaid traces, turn off the Show Detected Peaks command.	
Display Options > Show Legend	Displays a legend for the sample groups at the top of the view.	
Display Options > Legend Size	Specifies the number of legend lines that you want the application to display.	
	Selections: 0 to 10	
Display Options > Relative Intensity	Displays the <i>y</i> -axis scale as relative intensity (0 to 100%). The legend changes to Intensity [%].	
	The default <i>y</i> -axis scale is an absolute scale; the <i>y</i> -axis legend is intensity [counts].	
Display Options > Crop to Detected Peaks	Redraws only the chromatographic peaks of the displayed traces. Does not redraw the baseline portions of the traces as you select different table rows.	
Manual Peak Integration	Use to add a manual peak to a specialized trace. Adding a manual peak adds the Manual Peaks table to a result file (see "To manually integrate chromatographic peaks" on page 299).	
	Selecting a trace in the Specialized Traces table activates this command.	

Table 79. Shortcut menu commands for the Chromatograms view (Sheet 4 of 4)

Table 80 describes the traces that you can display in the Chromatograms view.

Trace type	Description	
UV	Displays a chromatogram created from the UV signal from a UV-Vis detector or the analog channel of a PDA detector.	
Analog	Displays a trace of response versus time.	
	Raw data files can contain analog data from a device that is hard-wired to the analog channels of a Thermo Scientific mass spectrometer.	
PDA total scan	Displays a chromatogram of the total absorbance for the entire scan wavelength range for each time point.	
PDA spectrum maximum	Displays a chromatogram of the highest absorbance reading in the wavelength range for each time point.	
PDA wavelength range	Displays a chromatogram of the total absorbance for the specified wavelength range for each time point.	
Base peak chromatogram (BPC)	Displays a chromatogram of the most intense mass spectral peak in the specified mass range for each time point.	
Total ion chromatogram (TIC)	Displays a chromatogram of the total intensity from all the mass spectral peaks in the specified mass range for each time point.	
Extracted ion chromatogram (XIC)	Displays an XIC trace, which is a mass range trace, when you select a row in any of these tables: Merged Features, Expected Compounds, Expected Compounds per File, Expected Compound Features, Compounds, Compounds per File, Unknown Compound Features, FISh Trace Fragments, or Specialized Traces.	
	The XIC trace is made up of the mass spectral peaks that match the specified mass value within the specified mass tolerance.	
Pattern trace	Displays a TIC trace of the summed intensities of the mass spectral peaks that match a specified pattern for each time point.	
FISh trace	Displays a TIC trace of the summed intensities of the mass spectral peaks in a fragmentation scan (MS/MS or MS ³) that match the predicted fragments of the selected library compound and its transformation products for each time point.	

Table 80. Chromatograms view traces

Working with the Mass Spectrum View

The Mass Spectrum view displays the spectral tree of a selected component in the result table and the mass spectrum of the selected scan in the spectral tree.

See these topics:

- Displaying a Mass Spectrum
- Changing the Zoom Level of the Mass Spectrum View
- Viewing Fragment Structures for Expected Compounds
- Creating a Mirror Plot
- Searching the mzCloud Library for a Matching Fragmentation Spectrum
- Spectral Tree Pane of the Mass Spectrum View
- Isotope Pattern Matching for Compounds with Formulas
- Mass Spectrum View Shortcut Menu Commands

Displaying a Mass Spectrum

To display a specific mass spectrum *

- 1. Open a result file.
- 2. If the Mass Spectrum view is closed, choose View > Mass Spectrum from the menu bar.
- 3. Do any of the following:
 - a. Select a row in any of these result tables.
 - Compounds Table

- Features Table
- Compounds per File Table
- Expected Features Table • mzCloud Results Table
- Expected Compounds Table
- Structure Proposals Table
- Expected Compounds per File Table The Mass Spectrum view displays the matching spectral tree and a zoomed-in view of

the full MS scan.

b. Select the scan of interest from the spectral tree.

Changing the Zoom Level of the Mass Spectrum View

In the Mass Spectrum view for a result file, you can independently zoom in or out on the x or y axis by dragging the cursor horizontally or vertically, respectively, or you can zoom in or out of a rectangular section of the plot.

Table 81. Working with the zoom level in the Mass Spectrum view

Task	Action
Zoom in on the <i>x</i> axis.	Drag the cursor to the right over the <i>m</i> / <i>z</i> range of interest.
Zoom out on the <i>x</i> axis.	Drag the cursor horizontally to the left over the m/z range of interest.
Zoom in on the <i>y</i> axis.	Drag the cursor vertically down the y axis.
Zoom out on the <i>y</i> axis.	Drag the cursor vertically up the <i>y</i> axis.
Zoom in on a section of the plot.	Drag the cursor diagonally across the section.
Undo the last zoom-in action.	Right-click the plot and choose Zoom Out .
Change the zoom level to the full m/z range on the x axis and the full response range on the y axis.	Right-click the plot and choose Undo All Zoom/Pan .

Viewing Fragment Structures for Expected Compounds

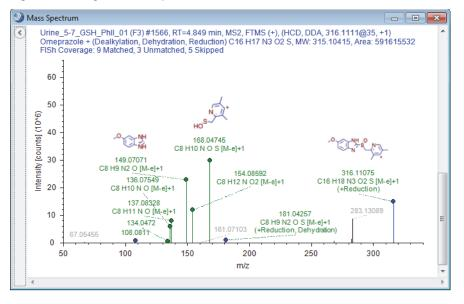
- To review the matching fragment structures predicted by the FISh Scoring node for expected compounds
- 1. Open a result file that was generated by a processing workflow that included the FISh Scoring node.
- 2. Select a compound in the Expected Compounds table.
- 3. (Optional) Enlarge the Mass Spectrum view.
- 4. In the spectral tree, select an MS2 or higher scan.
- 5. Review the color-coded mass spectral peaks, the theoretical fragment structures, and the transformations for the shifted mass spectral peaks.

The FISh Scoring Node annotates centroids that match the m/z value of a theoretical fragment ion with its theoretical structure and color-codes the centroids in a fragmentation scan as follows.

Color	Meaning
() Green	Direct match—Matches the m/z value of a theoretical fragment ion.
() Blue	Shifted match—Matches the m/z value of a theoretical fragment ion with at least one transformation applied.

Figure 87 shows an annotated fragmentation spectrum.





Creating a Mirror Plot

Create a mirror plot to compare an experimental spectrum to a reference spectrum.

To create a mirror plot of two fragmentation scans

- 1. Open a result file with fragmentation scans.
- 2. If the Mass Spectrum view is closed, choose View > Mass Spectrum from the menu bar.
- 3. In the spectral tree pane, select a reference scan.
- 4. Right-click the Mass Spectrum view and turn off the **Show Library Spectra as Reference** command if it is enabled.

Note The Use As Reference command is unavailable when the Show Library Spectra As Reference command is enabled.

5. Right-click the Mass Spectrum view and choose Use As Reference.

6. Select the fragmentation scan that you want to compare.

The reference scan appears on the bottom and the comparison scan appears on the top.

7. To remove the reference scan from the plot, right-click the Mass Spectrum view and choose **Clear Reference**.

Searching the mzCloud Library for a Matching Fragmentation Spectrum

* To manually search the mzCloud database for a matching fragmentation spectrum

Note The mzCloud database is only compatible with the Internet Explorer[™] web browser. To access the mzCloud spectral database from the Compound Discoverer application, set Internet Explorer as your default Internet browser.

- 1. Open a result file with data-dependent fragmentation data.
- 2. In the spectral tree pane of the Mass Spectrum view, select a fragmentation scan.
- 3. Right-click the spectrum plot and choose **Submit To mzCloud**.

The online mzCloud application opens to the Select Spectrum dialog box and displays the selected query spectrum.

4. Check the settings and click **OK**.

Spectral Tree Pane of the Mass Spectrum View

The collapsible pane on the left of the Mass Spectrum view contains a spectral tree with the high-resolution scans for preferred ions that elute within the following retention time window:

• Peak apex (RT) ± the peak's full width at half maximum (FWHM)

-or-

• From the start time to the end time for the chromatographic peak as defined by the peak detection algorithm

For MS1 scans, the tree lists the scan number, the retention time at the apex of the chromatographic peak, the mass analyzer, and the scan polarity.

For MSn scans, the tree lists the scan number, the retention time at the apex of the chromatographic peak, the mass analyzer, the scan polarity, and the fragmentation information, including the scan power, collision cell type, and the fragmentation type [data-dependent acquisition (DDA) or data-independent acquisition (DIA)].

Isotope Pattern Matching for Compounds with Formulas

Selecting a row in the following result tables populates the Mass Spectrum view with a spectral tree for the selected retention time in the left pane and the first MS1 scan for the selected retention time in the right pane.

- Expected Compounds
- Expected Compounds per File
- Expected Features
- Compounds

The MS1 scan shows the isotope pattern fit for the detected compound. Colored rectangles highlight the mass spectral peaks (centroids) that match the theoretical isotope pattern. These rectangles have a minimum display width to ensure that they are still visible when you zoom out or use the Undo All Zoom/Pan shortcut menu command.

Note The isotope pattern fit algorithm is "resolution aware"; that is, in addition to the list of elemental compositions provided by the Generate Expected Compounds node, it uses the resolution information provided with the scan data to perform an isotope pattern fit and calculate a spectral distance score.

If the resolution information is unavailable, it uses the setting for the Unrecognized MS Resolution parameter in the Select Spectra node.

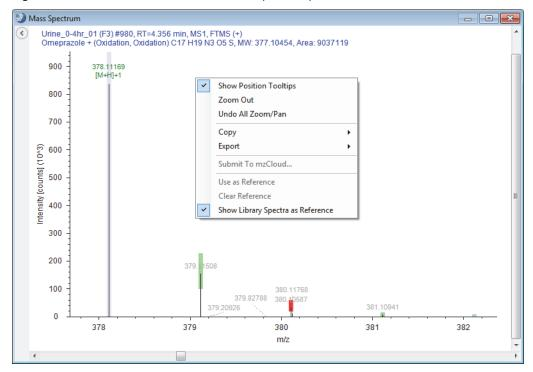


Figure 88. MS1 scan with color-coded mass spectrum peaks and the shortcut menu

Table 82 describes the color coding for the centroids in an MS1 spectrum.

Table 82. Color coding for the centroids in an MS1 spectrum for an expected compound (per file)

Color	Meaning
() Lavender	The labeled centroid matches the monoisotopic mass of the expected compound ion.
() Green	The labeled centroid matches the delta mass and the relative intensity of the theoretical isotope pattern within the specified tolerances.
() Red	The expected centroid for this m/z value is missing or its intensity does not fall within the tolerance range for the theoretical isotope pattern.
() Light blue	The expected centroid for this m/z value might be missing because its theoretical intensity is at the level of the baseline noise.

Mass Spectrum View Shortcut Menu Commands

Table 83 describes the shortcut (right-click) menu commands for the Mass Spectrum view (Figure 88 on page 309).

Command	Description	
Show Position Tooltips	Displays the <i>m</i> / <i>z</i> value and intensity of the mass spectrum peaks as you point to them.	
Zoom Out	Zooms out to the full scan range.	
Undo All Zoom/Pan	Zooms out to the full scan range.	
Copy > Image	Copies the mass spectrum as a bitmap (raster) image to the Clipboard.	
Copy > Points	Copies the data points and the scan header for the selected scan to the Clipboard. Also copies all text annotations, such as the FISh fragment annotations and the adduct information. Use this command to copy the FISh annotations to the Clipboard.	
Copy > Raw Points	Copies the data points for the selected scan to the Clipboard.	
	Use this command to copy points to a library search application.	
Export > Image As	Opens the Save As dialog box, where you can type a name and select a file type for the current scan. The available file types are BMP, EMT, GIF, JPG, PNG, and TIF.	
Export > Points As	Opens the Save As dialog box, where you can type a name and select a file type for the current scan. The available file type is TXT.	

Table 83. Mass Spectrum view shortcut menu commands (Sheet 1 of 2)

Command	Description
Submit to mzCloud	Opens the Spectrum Search dialog box for the mzCloud database where you can submit a search for the selected fragmentation scan.
	Available for an MS/MS data-dependent scan.
	IMPORTANT The mzCloud web application is only compatible with the Internet Explorer [™] web browser.
	Tip To test your computer's ability to connect to the mzCloud database, choose Help > Communication Tests from the application toolbar. Click the mzCloud tab, and then click Run Tests .
	To set up the mass tolerances for manual searches, choose Help > Configuration from the application toolbar. In the left pane, select Submit Single Spectrum to mzCloud Options . Edit the settings as applicable.
Use As Reference	Creates a mirror plot that initially consists of the currently selected scan. When you select another scan, the reference scan remains in the $-y$ -axis portion of the graph and the new scan appears in the $+y$ -axis portion of the graph.
	Unavailable when the Show Library Spectra Reference command is enabled.
Clear Reference	Removes the reference plot from the Mass Spectrum view.
Show Library Spectra as Reference	Displays the recalibrated library spectrum from the mzCloud database in the bottom portion of the mirror plot.
	Available when you select a row in the mzCloud Results table.

Table 83. Mass Spectrum view shortcut menu commands (Sheet 2 of 2)

Working with the Result Charts

Use the Result Charts view to plot the data in a result table as a scatter plot, histogram, bar chart, or pie chart.

Follow these topics as needed:

- Opening the Result Charts View
- Using the Copy, Export, and Zoom Commands for a Result Charts View
- Displaying and Pinning the Options Pane for the Result Charts View
- Plotting the Data as a Histogram Chart

- Plotting the Data as a Bar Chart
- Plotting the Data as a Pie Chart

Opening the Result Charts View

To open the Result Charts view

- 1. Open a result file.
- 2. In the menu bar, choose **View > Result Charts**.

By default, the Result Charts view opens as a floating window (Figure 89). You can resize the window, drag the window to another screen, or dock the window.

The Options pane for modifying the appearance of a chart view is a collapsible pane to the left of each chart.

Figure 89. Result Charts view with the collapsed (unpinned) Options pane at the left

🥑 F	🖉 Result Charts 📃 📼 💽		
Sca	atter Chart Histo	ogram Chart Bar Chart Pie Chart	
Options	Data Source: X Data:	ChemSpider Results Y Data:	Refresh Z Data:

Related Topics

- Opening, Closing, and Updating Result Files
- Rearranging the Tabbed Documents and Graphical Views

Using the Copy, Export, and Zoom Commands for a Result Charts View

* To access and use the shortcut menu commands for a Result Charts page

- 1. On the selected page of the Result Charts view, select the variables that you want to plot.
- 2. Click **Refresh** to plot the data.
- 3. Right-click the chart and choose one of the commands listed in Table 84.

Command	Availability	Function
Undo All Pan/Zoom	Scatter Chart, Histogram Chart, and Bar Chart	Undoes all panning and zooming. Returns the view to the default magnification.
Zoom Out	Scatter Chart, Histogram Chart, and Bar Chart	Undoes the last zoom-in action.
Copy > Image	All	Copies a raster image of the plot to the Clipboard.
Copy > Data	All	Copies a list of the data points to the Clipboard.
Export > Image As	All	Saves the plot in the selected file format.
		File types: EMF, PNG, GIF, JPG, TIF, and BMP
Export > Data As	All	Exports a list of data points to a text file.

Table 84. Common shortcut menu commands for the Result Chart views^a

^a The shortcut menu for the Scatter Chart view has additional commands.

Displaying and Pinning the Options Pane for the Result Charts View

The Options pane of the Result Charts view contains the formatting options for the grid lines, fonts, and so on.

To display and pin the Options pane

- 1. Point to the vertical Options tab on the left.
- 2. To keep the pane open, click the pin icon, +>, in the upper right corner of the Options pane.

Plotting the Data as a Histogram Chart

Use the Histogram Chart page of the Result Charts view to plot the frequency distribution of a variable (result table column). You can display the data as a column chart, a bar chart, a line, or a stepped line chart.

- To display the data as a histogram
- 1. Open a result file.
- 2. Choose View > Result Charts.

- 3. Click the Histograms Chart tab.
- 4. From the Data Source list, select the variable (result table column) that you want to plot.
- 5. Click Refresh.
- 6. (Optional) To change the appearance of the histogram, do the following:
 - a. Open the Options pane.
 - b. Modify the settings for the colors, labels, and legends of the display, as necessary.

-or-

Click **Load** to load the settings that you most recently saved.

c. To save the settings, click **Save** in the Options pane.

The new settings overwrite any previously saved settings. The application loads these new settings when you click Load.

d. To return the settings to the original default settings, in the Options pane, click **Factory Defaults**.

Table 85 describes the parameters for the Histogram Chart page.

Parameters	Description
Data Source	Specifies the source of the data that you want to plot. The available data sources are the result table columns in the result file.
Refresh	Refreshes the display with data points from the selected data source.
Options pane commands	
Load	Loads the Option pane settings that you saved with the Save command.
Save	Saves the settings that you selected in the Options pane.
Factory Defaults	Resets the settings of the options in the Options pane to the defaults in effect when you installed the application.

Table 85. Histogram Chart parameters (Sheet 1 of 6)

Parameters	Description
1. Chart Options	
Chart Type	Determines how the application plots the data.
	• (Default) Column: Displays the data as columns extending from bottom to top.
	• Bar: Displays the data as bars extending from left to right.
	• Line: Displays the data as a line. Reported and excluded items are not stacked as they are in column and bar types.
	• StepLine: Displays the data as a line drawn in a series of 90-degree angles.
Show Cumulative	Determines whether the application displays a cumulative histogram, which can be useful if you want to compare distribution curves. For these charts, the application calculates the column height as the height of the current column plus the sum of all previous columns. The rightmost column height is therefore the total count or 100%, depending on the <i>y</i> -axis settings.
	• (Default) True: displays a cumulative histogram.
	• False: Displays a non-cumulative histogram.
Horizontal Grid Lines	Determines whether the histogram displays horizontal grid lines and specifies the style of these lines.
	• (Default) None: Displays no horizontal grid lines on the histogram.
	• Solid: Displays solid horizontal grid lines on the histogram.
	• Dotted: Displays dotted horizontal grid lines on the histogram.
Vertical Grid Lines	Determines whether the histogram displays vertical grid lines and specifies the style of these lines.
	• (Default) None: Displays no vertical grid lines on the histogram.
	• Solid: Displays solid vertical grid lines on the histogram.
	• Dotted: Displays dotted vertical grid lines on the histogram.

 Table 85.
 Histogram Chart parameters (Sheet 2 of 6)

Parameters	Description
2. Column Options	
Show Column Amount	Determines whether the chart displays the amount or count of items.
	• True: Displays the amount or count of items in the chart.
	• (Default) False: Does not the display the amount or count or items in the chart.
Show Percentages	Determines whether percentage values appear above the columns or to the right of the bars.
	• True: Displays percentage values above the columns or to the right of the bars.
	• (Default) False: Does not display percentage values above the columns or to the right of the bars.
Column Display	Specifies the appearance of the columns or bars.
	• (Default) Flat: Displays the columns or bars as flat rectangle
	• Cylinder: Displays the columns or bars as cylinders.
	• Embross: Displays the columns or bars as three-dimensional rectangles.
	• LightToDark: Displays the columns or bars as shaded rectangles.
Column Width	Specifies the relative width of the columns as a fraction between and 1.
	Default: 0.8
Column Label Font	Specifies the font of the column labels that appear on top of the bars in the histogram. These labels are visible if you set Show Column Amount to True.
	Default: 8-point Arial
3. Axis Options	
X-Axis Number Format	Specifies the notation of the numbers used for the <i>x</i> axis.
	• (Default) Decimal: Uses decimal notation.
	• Scientific: Uses scientific notation.

Table 85. Histogram Chart parameters (Sheet 3 of 6)

Parameters	Description
X-Axis Title	Specifies the label for the <i>x</i> axis.
	The default <i>x</i> -axis title is the category selected from the Data Source list.
Y-Axis Type	Specifies the axis type (scale) of the y axis.
	• (Default) Linear: Plots the data on a linear scale.
	• Log: Plots the data on a logarithmic scale.
	• Percent: Plots the data as a percentage of the number of items.
Y-Axis Title	Specifies the label for the <i>y</i> axis. The default <i>y</i> -axis title is Count.
Reduce Number of Axis Labels	Determines whether the application increases readability by reducing the maximum number of axis labels to 30. If the chart includes more than 30 values, it displays only every second or every third label.
	• (Default) True: Reduces the maximum number of axis labels to 30.
	• False: Does not reduce the maximum number of axis labels to 30.
Axis Title Font	Specifies the font used to denote the labels of the x and y axes.
	Default: 12-point Arial
Axis Scale Font	Specifies the font used to denote the scale of the x and y axes.
	Default: 10-point Arial
4. Binning Options	
Binning Method	Specifies the number of data groups to display or the width of a single data group.
	• (Default) Auto: Groups the data by estimating the number of columns needed to display the number of data items.
	• FixedWidth: Groups the data according to the Width value.
	• FixedNumber: Groups the data according to the Number of Categories value.
Number of Bins	Specifies the number of categories used to group the data.
	Default: 20
	Note For discrete numbers, the actual group number might be different.

Table 85. Histogram Chart parameters (Sheet 4 of 6)

Parameters	Description
Bin Width	Specifies the width of a single category used to group the data.
	Default: 1
	Note For discrete numbers, the actual group number might be different.
Use Full Series Value Range	Determines the range of data values that the application uses to compile the histogram.
	• (Default) True: Uses all data for the histogram.
	• False: Uses only the data between the values specified by the Minimum Value option and the Maximum Value option.
Minimum Value	Specifies the minimum value of the displayed data range. When you use this parameter and the Maximum Value parameter, set Use Full Data Range to False. Use these two parameters when you want to show only a subrange of the data in a histogram.
Maximum Value	Specifies the maximum value of the displayed data range. When you use this option and the Minimum Value option, set the Use Full Data Range parameter to False.
5. Legend Options	
Show Legend	Determines whether a legend appears and where it appears.
	• (Default) None: Does not display a legend.
	• Top: Displays a legend at the top of the histogram.
	• Left: Displays a legend to the left of the histogram.
	• Bottom: Displays a legend at the bottom of the histogram.
	• Right: Displays a legend to the right of the histogram.
Legend Font	Specifies the font for the legend.
Legend Font	-F

 Table 85.
 Histogram Chart parameters (Sheet 5 of 6)

Parameters	Description
6. Series Options	
Show Only Checked Items	Determines whether the chart is compiled from data in all result rows or only data in result rows marked by check marks.
	• True: Compiles the chart only from data in result rows marked by check marks.
	• (Default) False: Compiles the chart from data in all result rows.
Target Series Color	Specifies the color of the target series.
	Default: Firebrick

Table 85. Histogram Chart parameters (Sheet 6 of 6)

Related Topics

• Displaying and Pinning the Options Pane for the Result Charts View

Plotting the Data as a Bar Chart

Bar charts plot categorical and ordinal data types in columns with the count of the data types as the column height.

The Data Source list contains the numerical data categories that are available for the bar chart. The Options pane contains the different options that you can use to customize the bar chart. Moving the cursor over the columns in the chart activates a tooltip with information about the data category.

To display the data as a bar chart

- 1. In an open result file, choose View > Result Charts.
- 2. Click the Bar Charts tab in the Result Charts view.
- 3. From the Data Source list, select the type of data to display as a bar chart.
- 4. Click **Refresh** to draw the chart.
- 5. (Optional) To change the chart's appearance, do the following:
 - a. Open the Options pane and adjust the colors, labels, and legends of the display. Or, click **Load** to load the option settings that you most recently saved.
 - b. To save all the settings, click **Save** in the Options pane.

The new settings overwrite any previously saved settings. The application loads these new settings when you click Load.

c. To return the Options pane settings to the original default settings, in the Options pane, click **Factory Defaults**.

Table 86 describes the parameters on the Bar Chart page of the Result Charts view.

Table 86. Bar Chart parameters (Sheet 1 of 3)

Command or option	Description
Options pane	
Load	Loads the Option pane settings that you saved with the Save command.
Save	Saves the settings that you selected in the Options pane.
Factory Defaults	Resets the options in the Options pane to the default settings in effect when you installed the application.
Chart Options	
Chart Type	Determines how the application plots the data.
	• (Default) Column: Displays the data as columns extending from bottom to top.
	• Bar: Displays the data as bars extending from left to right.
Horizontal Grid Lines	Determines whether the bar chart displays horizontal grid lines and specifies the style of these lines.
	• (Default) None: Displays no horizontal grid lines in the bar chart.
	• Solid: Displays solid horizontal grid lines in the bar chart.
	• Dotted: Displays dotted horizontal grid lines in the bar chart.
Axis Options	
X-Axis Title	Specifies the label for the <i>x</i> axis. The default <i>x</i> -axis title is the category selected from the Data Source list.
Y-Axis Title	Specifies the label for the y axis. The default y -axis title is Count.
Y-Axis Type	Specifies the axis type (scale) of the y axis.
	• (Default) Linear: Plots the data on a linear scale.
	• Log: Plots the data on a logarithmic scale.
	• Percent: Plots the data as a percentage of the number of items.
Axis Title Font	Specifies the font used to denote the labels of the <i>x</i> and <i>y</i> axes.
	Default: 12-point Arial

Command or option	Description
Axis Scale Font	Specifies the font used to denote the scale of the x and y axes.
	Default: 10-point Arial
Bar Options	
Show Rotated Scale	Determines whether the scale labels along the <i>x</i> axis are slightly rotated to the right.
	• (Default) True: Rotates the scale labels.
	• False: Does not rotate the scale labels.
X-Axis Scale Angle	Specifies the angle between the label and the <i>x</i> axis for the labels when the Show Rotated Labels parameter is set to True.
	Range: –90 to 90 degrees
	Default: 30
Bar Display	Specifies the appearance of the columns or bars.
	• (Default) Flat: Displays the columns or bars as flat rectangles
	• Cylinder: Displays the columns or bars as cylinders.
	• Embross: Displays the columns or bars as three-dimensional rectangles.
	 LightToDark: Displays the columns or bars as shaded rectangles.
Bar Width	Specifies the relative width of the columns.
	Range: 0.1–1.0
	Default: 0.8
Show Amount	Determines whether the amount, or count of items, is displayed in the chart.
	• True: Displays the amount or count of items in the chart.
	• (Default) False: Does not the display the amount or count of items in the chart.

 Table 86.
 Bar Chart parameters (Sheet 2 of 3)

Command or option	Description
Show Percentage	Determines whether percentage values are displayed above the columns or to the right of the bars.
	• True: Displays percentage values above the columns or to the right of the bars.
	• (Default) False: Does not display percentage values above the columns or to the right of the bars.
Column Label	Specifies the font of the column labels that appear on top of the bars in the bar chart.
	Default: 8-point Arial
Legend Options	
Show Legend	Determines whether a legend appears and where it appears.
	• (Default) None: Does not display a legend.
	• Top: Displays a legend at the top of the bar chart.
	• Left: Displays a legend to the left of the bar chart.
	• Bottom: Displays a legend at the bottom of the bar chart.
	• Right: Displays a legend to the right of the bar chart.
Legend Font	Specifies the legend font.
	Default: 8-point Arial
Series Options	
Show Only Checked Items	Determines whether the chart is compiled from data in all result rows or only data in result rows marked by check marks.
	• True: Compiles the chart only from data in result rows marked by check marks.
	• (Default) False: Compiles the chart from data in all result rows.
Series Color	Specifies the color of the bars.
	Default: CornflowerBlue
Data Source	Displays the result category used to plot the data.

Table 86. Bar Chart parameters (Sheet 3 of 3)

Plotting the Data as a Pie Chart

The Pie Chart page shows several categories of data as a solid circle composed of slices (a pie) or as a ring (a doughnut). You can use a pie chart or a doughnut chart to indicate the relative size of quantities of data.

✤ To display the data as a pie chart

- 1. With an active result file, choose **View > Result Charts** from the menu bar.
- 2. Click the **Pie Chart** tab in the Result Charts view.
- 3. From the Data Source list, select the type of data to display.
- 4. Click **Refresh** to draw the chart.
- 5. (Optional) To change the chart's appearance, do the following:
 - a. Open the Options pane and adjust the colors, labels, and legends of the display and specify how the chart displays small slices. Or, click **Load** to load the option settings that you most recently saved.

Tip By default, the application consolidates small slices of 5% or less. To change this setting, do one of the following:

• Select False for Collect Small Segments.

-or-

- Change the Small Slice Threshold (%) setting.
- b. To save all the settings, click **Save** in the Options pane.

The new settings overwrite any previously saved settings. The application loads these new settings when you click Load.

When you close the Result Charts view, the application stores the chart settings that you selected. When you reopen the chart, it displays these stored settings if they are available.

c. To return the Options pane settings to the original default settings, in the Options pane, click **Factory Defaults**.

Table 87 describes the parameters on the Pie Chart page of the Result Charts view.

Table 87. Pie Chart page parameters (Sheet 1 of 4)

Command or Option	Description
Options pane	
Load	Loads the Option pane settings that you saved with the Save command.
Save	Saves the settings that you selected in the Options pane.

Command or Option	Description
Factory Defaults	Resets the options in the Options pane to the default settings in effect when you installed the application.
Chart Options	
Chart Type	Specifies the type of chart to display:
	• (Default) Pie: Displays the chart as a solid circle composed of slices.
	• Doughnut: Displays the chart as a ring.
Vertical Rotation	Specifies the angle of rotation around the vertical axis.
	Default: 0
Angle of First Slice	Specifies the location of the first chart slice.
	Default: 0
Pie Slice Options	
Show Type Name with Value	Determines whether the application displays the name of the data group next to the value.
	• True: Displays the name of the data group next to the value.
	• (Default) False: Does not display the name of the data group next to the value.
Show Slice Amount	Determines whether the application displays the amount of each slice.
	• (Default) True: Displays the amount of each slice.
	• False: Does not display the amount of each slice.
Show Slice Percentage	Determines whether the application displays the percentage of each slice.
	• (Default) True: Displays the percentage of each slice.
	• False: Does not display the percentage of each slice.
Label Style	Specifies the label style of the chart segments.
	• Disabled: Does not display a label.
	• (Default) Inside: Displays the label on top of the chart segment.
	• Outside: Displays the label outside the chart segment.

Table 87. Pie Chart page parameters (Sheet 2 of 4)

Command or Option	Description
Labels Font	Specifies the font of the data set labels.
	Default: 9-point Microsoft Trebuchet bold
Small Slices Options	
Collect Small Segments	Determines whether the application consolidates small segments together into a single slice.
	• (Default) True: Consolidates small segments into a single slice.
	• False: Leaves small segments as is.
Small Slice Threshold (%)	Specifies a percentage threshold for the slices to be consolidated into a single slice.
	Default: 5%
Small Slice Collection Label	Specifies the label for the pie segment composed of consolidated segments.
	Default: Other
Small Slice Collection Color	Specifies the color of the pie segment composed of consolidated segments.
	Default: Gray
Show as Supplemental Pie	Determines whether to display small segments as a supplemental pie chart.
	• True: Displays small segments as a supplemental pie chart.
	• (Default) False: Leaves small segments as is.

Table 87. Pie Chart page parameters (Sheet 3 of 4)

Command or Option	Description
Supplemental Pie Size	Specifies the size of the supplemental pie chart relative to the original pie chart.
	• Largest: Displays the supplemental pie chart as much larger than the original pie chart.
	• Larger: Displays the supplemental pie chart as larger than the original pie chart.
	• Comparable: Displays the supplemental pie chart as about the same size as the original pie chart.
	• (Default) Smaller: Displays the supplemental pie chart as smaller than the original pie chart.
	• Smallest: Displays the supplemental pie chart as much smaller than the original pie chart.
Data Source	Displays the result category used to plot the data.

Table 87. Pie Chart page parameters (Sheet 4 of 4)

Working with a Scatter Plot

Use the Scatter Chart page to set up a scatter plot for visualizing whether there is a linear or logarithmic relationship between two or three variables (columns) in a result table.

To plot the data as a scatter chart, follow these topics as needed:

- Setting Up a Scatter Plot
- Using a Filter Set to Filter the Scatter Plot
- Customizing the Appearance of a Scatter Plot
- Customization Options for a Scatter Plot
- Scatter Chart Parameters
- Scatter Chart Shortcut Menu Commands

Related Topics

• Working with the Result Charts

Setting Up a Scatter Plot

✤ To set up a scatter plot

- 1. Open a result file.
- 2. In the menu bar, choose **View > Result Charts**.

By default, the Result Charts view opens as a floating window.

- 3. Click the **Scatter Chart** tab.
- 4. In the Data Source list, select one of the available result tables.

The available selections depend on the workflow nodes in the processing workflow.

- 5. Select the variables as follows:
 - a. In the X Data list, select the variable that you want to plot against the *x* axis.
 - b. In the Y Data list, select the variable that you want to plot against the *y* axis.
 - c. To create a three-dimensional scatter plot, select the variable in the Z Data list that you want to plot against the *z* axis.

Selecting a data value for the z axis adds a color gradient to the plotted data points ranging from the lowest to the highest Z data value.

6. Click Refresh.

Depending on your selections, a two- or three-dimensional linear plot of the data points appears. With the default appearance settings, the data points appear as blue circles in a 2D plot and as circles of varying colors in a 3D plot. For a 3D plot, a color legend for the lowest to the highest Z data value appears to the right of the scatter chart (Figure 90).

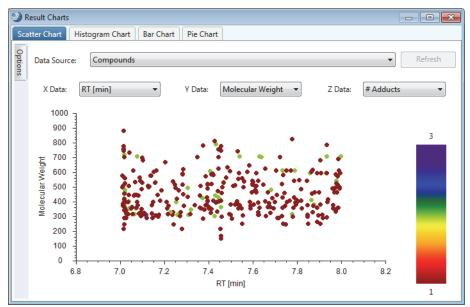


Figure 90. Three-dimensional scatter chart

Using a Filter Set to Filter the Scatter Plot

- * To interactively filter the scatter plot by using the Result Filters view
- 1. Set up and apply a set of result filters in the Result Filters view.

In the Scatter Chart view, the Refresh button turns orange.

2. Click **Refresh** to refresh the scatter chart plot.

Customizing the Appearance of a Scatter Plot

In the Options pane, do any of the following:

• To change the scaling, colors, labels, and legends in the display, under Axis Options, make the appropriate changes.

The application applies the changes as you make them.

• To change the font size or font type of the axis labels, click the expand icon to the left of Axis Scale Font or Axis Title Font. Then, make the appropriate selections or click the browse icon to open the Font dialog box where you can make your selections.

A browse icon and a set of font parameters appear.

• To return the option settings to the original default settings, click Factory Defaults.

Related Topics

• Displaying and Pinning the Options Pane for the Result Charts View

Customization Options for a Scatter Plot

Table 88 describes the formating options for the Scatter Chart page.

Table 88. Options pane for a Scatter Chart

Parameter	Description
Buttons	
Load	Loads the saved Options pane settings. Only click Load to apply a set of saved settings; otherwise, the view reverts to the last saved set.
Save	Saves the new settings.
Options pane. Howe	rt view automatically updates as you change the settings in the ver, the application does not save the settings until you click Save. If e saving the new settings, the view reverts to the previously saved
Factory Defaults	Resets the options in the Options pane to the default settings in effect when you installed the Compound Discoverer application.
Axis Options	
X Axis Type	Specifies the axis type (scale) of the <i>x</i> axis: Linear or Logarithmic.
Y Axis Type	Specifies the axis type (scale) of the <i>y</i> axis: Linear or Logarithmic.
Z Axis Type	Specifies the axis type (scale) of the <i>z</i> axis: Linear or Logarithmic.
Axis Scale Font	Specifies the font used to denote the scale of the x and y axes.
Axis Label Font	Specifies the font used for the titles of the <i>x</i> and <i>y</i> axes.
X-Axis Title	Specifies the title of the <i>x</i> axis.
Y-Axis Title	Specifies the title of the y axis.
Series Options	
Points	Specifies the appearance of the points in the scatter chart. Also specifies whether to show the points in the scatter chart.
Selected Points	Specifies both the appearance of the selected points in the scatter chart and whether to show the points in the scatter chart.
Filtered-out Points	Specifies the appearance of the filtered-out points in the scatter chart and whether to show the points in the scatter chart.
Excluded Points	The Compound Discoverer workflow nodes do not generate excluded data points.

Scatter Chart Parameters

Table 89 describes parameters that are visible on the Scatter Chart page. Use these parameters to set up the scatter plot.

Table 89. Scatter Chart parameters

Parameter	Description	
Data Source	Specifies the data source for the plot. The data source is one of the available result tables produced by the processing workflow.	
Note The X, Y, and Z Data boxes list the available variables for the selected data source. The variables are the available columns in the selected data source (result table).		
X Data	Specifies the variable to plot against the <i>x</i> axis.	
Y Data	Specifies the variable to plot against the <i>y</i> axis.	
Z Data	Specifies the variable to plot against the <i>z</i> axis.	
Plot grid	Two-dimensional grid where the application plots the data points. By default, the plot area has no grid lines.	
	To add horizontal and vertical lines to the plot, make the appropriate selection under Options in the Options pane.	
Axis labels	The default axis labels are the selected variable names.	
Color legend	When you create a 3D plot, the scatter chart includes a color legend for the <i>z</i> -axis color gradient. The numeric value of the highest <i>z</i> -axis data point appears above the color legend.	
Options panel	Use the parameters in this pane to customize the Scatter Chart page. For more information, see Table 88.	
Buttons		
Refresh	Refreshes the content of the chart area.	

Scatter Chart Shortcut Menu Commands

Command	Description
Show Position Tooltips	Displays a tooltip when you place the cursor over a data point.
Zoom Out	Decreases the zoom of both axes.
Undo All Zoom/Pan	Displays the full <i>x</i> -axis range of the plot.
Copy > Image	Copies the scatter chart plot as a bitmap (raster) image to the Clipboard.
Copy > Data	Copies the data points for the selected variables to the Clipboard.

Table 90. Scatter Chart shortcut menu commands (Sheet 1 of 2)

Command	Description
Export > Image As	Saves the plot to any of these file types: EMF, PNG, GIF JPG, TIF, and BMP
Export > Data As	Opens the Save As dialog box where you can save the data as a plain text file.
	By default, the application saves the file to the last open folder and uses the following convention to name the file:
	ResultTable_x_XDataSelection_Range_YDataSelection_Range _z_ZDataSelection_Range_DataPointType.txt
	Use the Browse icon to select a different folder.
	Exporting data from a scatter plot creates three text files: <i>File name</i> Filtered-out Points.txt, <i>File name</i> Points.txt, and <i>File name</i> Selected Points.txt.
Select Item for Point	Highlights the appropriate row in the result table (selected Data Source).
Check Point	By default, changes the selected point to a red diamond and selects the check box in the Checked column for the selected point (row in the selected Data Source result table). You can change the appearance of selected points by making the appropriate selections in the Series Options > Checked Points area of the Options pane. When you save the data to a text file, the Selected Points.txt file lists the selected points.
	To select a point, right-click the point of interest on the plot and choose Check Point .
Uncheck Point	Undoes the selection of a selected point. You can change the appearance of points by making the appropriate selections in the Selected Options > Points area of the Options pane.
Check All Visible Points	Selects all of the visible points in the scatter plot.
Uncheck All Visible Points	Undoes the selection of the visible points in the scatter plot.

Table 90. Scatter Chart shortcut menu commands (Sheet 2 of 2)

Working with the Trend Chart View

Use the Trend Chart view to compare the chromatographic peak areas for compounds by sample group.

You define the sample groups by selecting one or more check boxes under Group By in the collapsible pane to the left of the chart. In addition, you can change the sort order of each sample group and the hierarchy of the sample groups.

You can use the Trend Chart view to plot the data for a compound (or feature) in these result tables: Compounds, Expected Compounds, and Merged Features. For a single compound, the chart can plot the data as a trendline plot or a box-and-whisker plot. For two or more compounds, the chart displays the data only as a trendline plot, with one trendline for each compound.

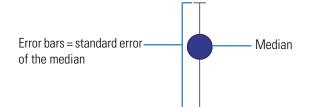
Note The trendline plot can plot sample groups that include only one data point; however, to plot the error bars for a group, it requires a minimum of two data points. With two data points, the circle represents the calculated median and the error bars represent the minimum and maximum areas.

The box-and-whisker plot requires a minimum of two data points to plot the box for a sample group. It does not plot sample groups that include only one data point. If none of the sample groups includes the minimum number of data points, the following text appears in the chart area: No results available to plot.

By default, the trendline plot type plots the median peak area for the selected component on the *y* axis against equally spaced sample groups on the *x* axis and connects the data points with straight lines. Each data point appears as a solid circle with two error bars. The circle represents the median (Figure 91). If the group includes a least three data points, the application uses the following equation to calculate the standard error of the median and draw the error bars:

standard error of the median = $\sqrt{\pi/2} \times \text{std. dev}(x)/(\sqrt{N})$

Figure 91. Trendline display of the data distribution

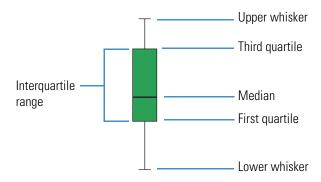


The box-and-whisker chart plots the peak area for a selected component on the y axis as a rectangle against equally spaced sample groups on the x axis. The height of the rectangle represents the peak areas in the interquartile range (Figure 92). The application uses the following equations to calculate the upper and lower whiskers:

Interquartile range (IQR) = Quartile 3 (Q3) – Quartile 1 (Q1) Upper whisker = Q3 + IQR × 1.5 Lower whisker = Q1 – IQR × 1.5

When the data set contains a small number of data points, the whiskers typically end at the highest and lowest data points. If the data set does not include a data point between the top of the interquartile range and the calculated value for the upper whisker, the application does not draw an upper whisker. If the data set does not include a data point between the bottom of the interquartile range and the calculated value for the lower whisker, the application does not draw a lower whisker.

Figure 92. Box-and-whisker display of the data distribution



Note To calculate the quartiles, the application uses a method that is similar to the type 6 method in the R statistical computing software.

Follow these topics as needed:

- Opening the Trend Chart View
- Defining the Sample Groups to Compare
- Comparing the Peak Areas for a Single Compound by Sample Group
- Changing the Sort Order of the Defined Groups
- Comparing the Peak Areas for Multiple Compounds by Group
- Changing the Hierarchy of the Variables Used for Grouping
- Showing the Error Bars in a Trendline Chart
- Trend Chart View Parameters

Opening the Trend Chart View

To open the Trend Chart view

- 1. Open the result file of interest.
- 2. From the menu bar, choose View > Trend Chart.

The Trend Chart view opens as a docked window to the right of the result tables.

- If the active result table does not contain a consolidated compounds list, the following text appears in the graph area: No Results Available to Plot. The Compounds, Expected Compounds, and Merged Features tables contain a consolidated compounds list.
- If the active result table contains a consolidated compounds list and you select a row, a box-and-whisker plot appears in the graph area with data from the first table row.

Defining the Sample Groups to Compare

- To define the sample groups for a trend chart plot
- 1. Open the result file of interest, and then open the Trend Chart view.
- 2. In the left pane, under Group By, select the appropriate check boxes to define the sample groups.

Comparing the Peak Areas for a Single Compound by Sample Group

- $\boldsymbol{\ast}$ To compare the peak areas for a single compound by sample group
- 1. Open a result file, open the Trend Chart view, and define the sample groups that you want to compare.
- 2. In the Compounds, Expected Compounds, or Merged Features table, select the compound (or feature) of interest.
- 3. Right-click the graph area and choose **Show Legend**.

The legend displays the sample group colors.

4. In the Plot Type list, select Trendline Chart or Box Whisker Chart.

Depending on the selection, either a trendline plot or a box-and-whisker plot appears in the graph area. A tooltip opens when you place the cross-hair cursor anywhere on a box or whisker in the box-and-whisker plot or on data point in a trendline plot.

Figure 93 shows a box-and-whisker plot. Placing the cross-hair cursor on a box or whisker opens a tooltip with descriptive statistics.

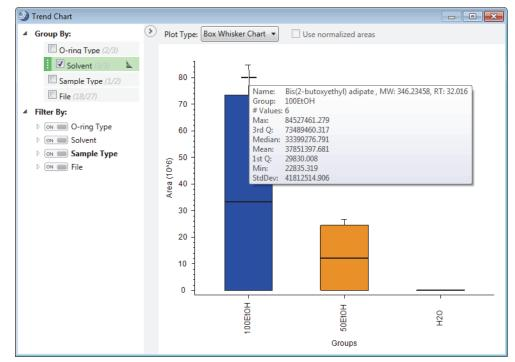


Figure 93. Box-and-whisker plot for one compound

Changing the Sort Order of the Defined Groups

In the Trend Chart view, you can display the defined groups from left to right in ascending or descending order by chromatographic peak area.

* To change the sort order of the defined groups in the Trend Chart view

In the left pane, under Group By, click the sorting icon, \blacksquare , next to the study variable to sort and choose **Sort Ascending** or **Sort Descending**.

Changing the sort order changes the order of the groups on the *x* axis (Figure 94).

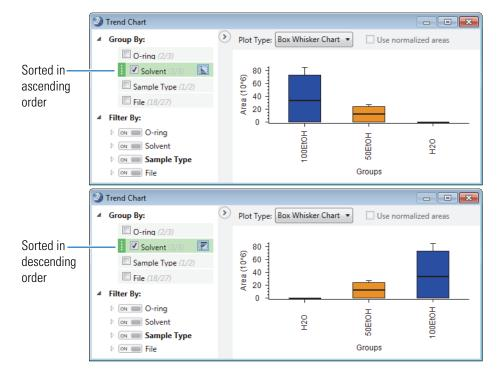


Figure 94. Trend chart with compounds grouped by solvent and sorted in ascending or descending order by chromatographic peak area

Comparing the Peak Areas for Multiple Compounds by Group

* To compare the peak areas for multiple compounds by group

- 1. Open the result file of interest, open the Trend Chart view, and define the sample groups.
- 2. To select the compounds to plot, press the CTRL key and select rows in the result table (Compounds, Expected Compounds, or Merged Features), taking care to avoid clicking an editable table cell.

The plot changes to a scaled trendline. The data points represent the group median. The legend displays the name (if available), molecular weight, and retention time of each selected compound.

Figure 95 shows a scaled trendline chart. To view descriptive statistics for the data points, click anywhere in the plot to activate the cross-hair cursor, and then place the cross-hair cursor on each data point of interest.

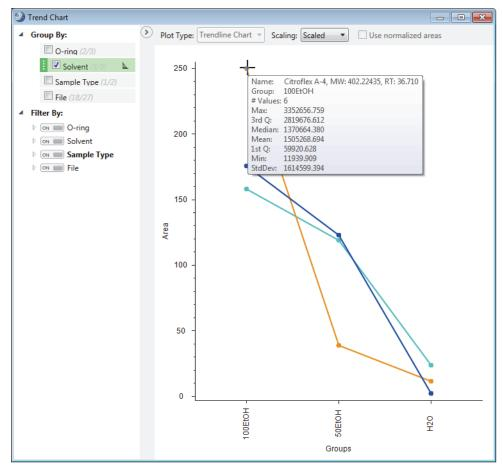


Figure 95. Trendline chart with three compounds

Changing the Hierarchy of the Variables Used for Grouping

Changing the hierarchy of the variables used for grouping in the Trend Chart view modifies the sample grouping.

* To change the hierarchy of the variables used for grouping

Use the handle () next to the variable to drag the variable up or down in the list.

Showing the Error Bars in a Trendline Chart

- ✤ To view error bars for each data point in a trendline chart
- 1. In the Scaling list above a trendline plot in the Trend Chart view, select Unscaled.
- 2. Right-click the plot and choose Show Standard Errors.

The application plots the group median with error bars for the standard error of the median (Figure 96).

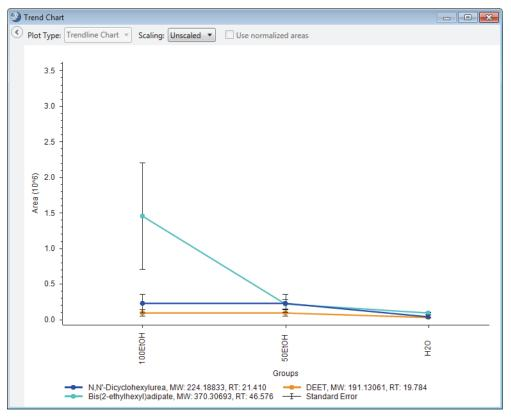


Figure 96. Trendline chart with error bars

Trend Chart View Parameters

Table 91 describes the parameters in the Trend Chart view.

Table 91. Trend Chart view parameters

Parameter	Description			
Plot Type	Controls the plot type in the graph area.			
	Selections: Trendline Chart or Box Whisker Chart			
	The Box Whisker Chart selection is only available for displaying a single compound.			
Scaling (for the trendline chart)	 Select the chart scaling: Unscaled—Enables the Show Standard Errors command in the shortcut menu. Scaled Scaled to Study Factor Value 			
Use Normalized Areas	Displays the normalized areas for the data.			
	Available when the processing workflow included the Normalize Areas node.			

Working with the Mass Defect Plot View

Use the Mass Defect Plot view to find similar compounds by visualizing the calculated mass defects for a set of compounds against their molecular weights. Your choices for calculating the mass defect in this view are the same as those for the Calculate Mass Defect and Filter By Mass Defect nodes.

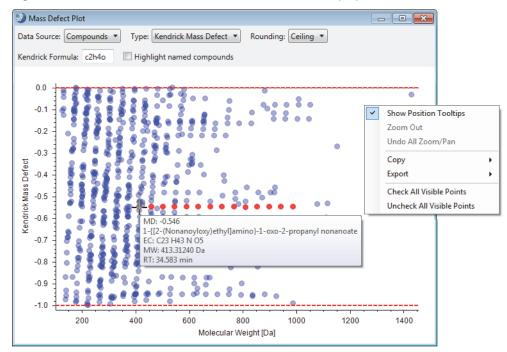
- * To set up the mass defect plot and examine it for similar compounds
- 1. Open a result file from a processing workflow that included the Detect Compounds node.
- 2. From the application menu bar, choose View > Mass Defect Plot.

The compounds listed in the Compounds table appear as blue circles in the Mass Defect plot. The *y*-axis label displays the mass defect type, and the red dashed lines indicate the valid range of the mass defect values.

- 3. To specify how to calculate the mass defect, do the following:
 - From the Type list, select the mass defect calculation. Then, if you selected Kendrick Mass Defect, enter the Kendrick formula.
 - From the Rounding list, select Ceiling, Floor, or Round.

- 4. To highlight the compounds that have an assigned name in orange, select the **Highlight** Named Compounds check box.
- 5. To work interactively with the plot, do the following:
 - Use the view's shortcut menu commands to do any of the following:
 - Zoom in or out of the plot.
 - Copy the image or the data points to the Clipboard.
 - Export the image or the data points to an external file.
 - Check or uncheck the visible data points. Checking a data point highlights the data point in red and places a check mark in the Checked column for the compound in the Compounds table.
 - Double-click a data point to navigate to the compound in the Compounds table.
 - Point to a compound to display a tooltip with information about the compound's mass defect, name, elemental composition, molecular weight, and retention time.

Figure 97. Mass Defect Plot view with its shortcut menu displayed



Related Topic

• Calculating the Mass Defect of an Elemental Composition

Working with the Principal Component Analysis View

Use the Principal Component Analysis view to visualize the correlation between multivariate data in a set of observations. A principal component analysis transforms a set of observations for possibly correlated variables into an artificial set of independent linear combinations of the original variables known as principal components (PC1, PC2, PC3 and so on). PC1 has the most variation and the highest principal component has the least variation.

The Principal Component Analysis view contains three pages: Scores Plot, Loadings Plot, and Variances Plot. The scores plot shows the correlation among the observations. The loadings plot shows the relationship among the variables for a given pair of principal components. The variance plot shows the percentage and cumulated percentage of the variance that a principal component accounts for. In general, as the proportion of variance increases for the first two or three principal components, the dissimilarity between the sample groups increases.

To work with the Principal Component Analysis view, follow these topics:

- Setting Up a Principal Component Analysis
- Interpreting the Scores Plot
- Interpreting the Loadings Plot
- Working Interactively with the Loadings Plot
- Interpreting the Variance Plot
- Principal Component Analysis Parameters

Setting Up a Principal Component Analysis

To set up a principal component analysis

- 1. Open a result file that contains any of these result tables: Compounds, Expected Compounds, or Merged Features.
- 2. From the menu bar, choose View > Principal Component Analysis.

If the result file contains a Compounds table, the Principal Component Analysis view opens with the Compounds table selected as the data source.

- 1. In the Data Source list, select the **Compounds**, **Expected Compounds**, or **Merged Features** table as appropriate for the data set.
- 2. In the X Data list, select PC1, PC2, PC3, PC4, or PC5.

In most cases, select PC1 or PC2, as these principal components have the most variation.

3. If the data points differ by several orders of magnitude, select the **Center and Scale** check box.

- 4. If the processing workflow included the Normalize Areas node, select the **Use Normalized Areas** check box as appropriate.
- 5. To display a legend at the bottom of the plot, right-click the plot and choose **Show** Legend.
- 6. To show a tooltip when you place the cross-hair cursor on a data point in the plots, right-click the plot and choose **Show Position Tooltips**.
- 7. To remove points from the plot, clear the check boxes under Filter By as appropriate.

Using the new population, the application recalculates the principal components, including their contribution to the variance, and shifts the coordinates of the remaining data points.

Interpreting the Scores Plot

Use the scores plot of the Principal Component Analysis view to interpret the relationship among the sample groups. Sample groups that are near each other are similar.

The Principal Component Analysis view is available for result files that include any of these tables: Compounds, Expected Compounds, or Merged Features.

To review and interpret the scores plot

- Check the percentage values for the principal components. The labels on the *x* and *y* axes include the proportion of variance that the principal components add to the total variance as a percentage.
- Place the cross-hair cursor on a data point to view its coordinates.

Figure 98 shows the scores plot for the compounds that leach out of four o-ring types soaked in ethanol. The black, brown, and white o-rings show a similar variance for PC1, whereas the red o-rings show a significant variance in the other direction for PC1. The brown and white o-rings show a similar variance for PC2.

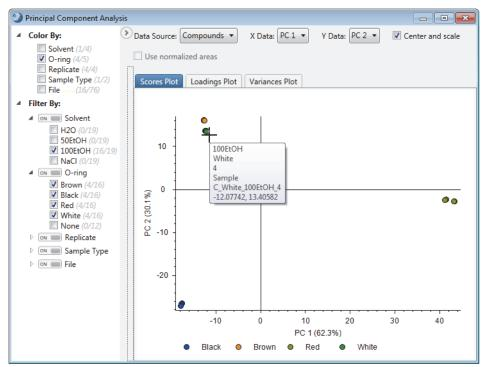


Figure 98. Scores plot for four o-ring types (white, brown, red, and black)

Interpreting the Loadings Plot

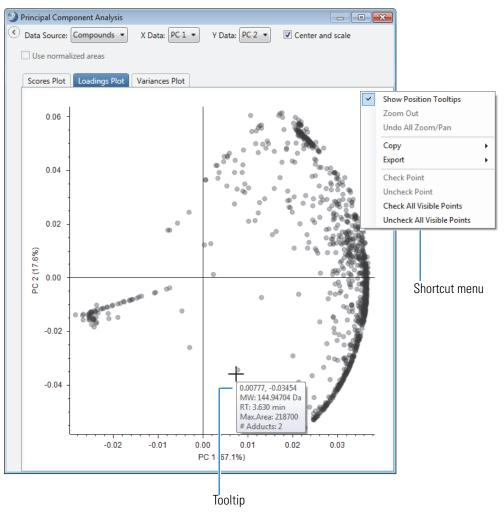
Use the loadings plot to interpret the relationship among the variables.

✤ To interpret the loadings plot

- 1. To determine the relative correlation between various data points, review their relative location in the plot.
 - Data points that are near each other are similar.
 - Data points that are on opposite sides of the origin have a negative correlation.
 - Data points in the corners of the plot have a strong contribution to both principal components—that is, these data points differentiate between groups.

2. Place the cross-hair cursor on a data point to display a tooltip with the following information about the compound: the principal component coordinates, molecular weight, retention time, maximum peak area, and number of adduct ions (Figure 99).

Figure 99. Loadings plot with its shortcut menu displayed



Working Interactively with the Loadings Plot

The Loadings Plot is interactive with the selected data source table—that is, double-clicking a data point in the loadings plot selects its row in the data source table and updates the other opened graphical views. In addition, checking data points (compounds, expected compounds, or features) in the Loadings Plot checks the corresponding data points in these interactive views—Differential Analysis, Partial Least Squares Discriminant Analysis, and Descriptive Statistics.

- * To select a data point or check one or more data points in the Loadings Plot
 - To select a data point, double-click it.

The application highlights the corresponding row in the data source table and updates the other opened and interactive views.

- To check a single data point, right-click the point and choose Check Point.
- To check multiple points, drag the cursor across a rectangular area of the plot to zoom in on that area, then right-click the plot and choose **Check All Visible Points**.

The application selects the check boxes for the checked data points in the data source table and changes the color of the selected points to blue in the Loadings Plot (Figure 100). The color of the checked data points also changes to blue in the interactive views.

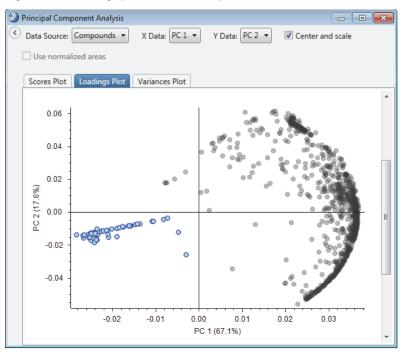


Figure 100. Loadings plot with checked points shown in blue

• To clear all the selected check boxes in the result table, right-click the result table and choose **Uncheck All > In This Table**.

Interpreting the Variance Plot

Use the variance plot to determine the relative differentiation between sample groups.

To interpret the variance plot

Compare the contribution of the first two or three principal components to the cumulative variance.

For example, Figure 101 shows the variance plot for the compounds leached from red o-rings soaked either in water or an aqueous solution of sodium chloride, and Figure 102 shows the variance plot for the same type of o-rings soaked either in water or ethanol.

The PC1–PC3 components make up more of the cumulative variance when comparing the extraction strength of water to ethanol than when comparing water to a salt solution. From these results, you can infer that the extraction strengths of water and ethanol differ more than the extraction strengths of water and a salt solution.

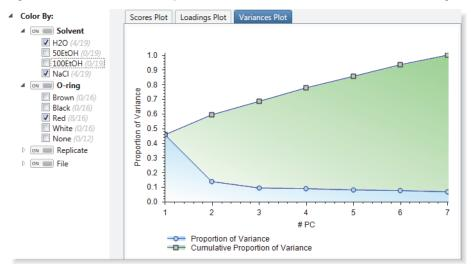
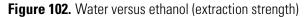
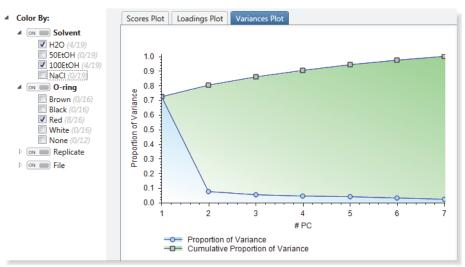


Figure 101. Water versus an aqueous solution of sodium chloride (extraction strength)





Principal Component Analysis Parameters

Table 92 describes the parameters in the Principal Component Analysis view.

-			
Parameter	Description		
Data Source	Specifies the source of the data. The selection list depends on the processing workflow.		
	Selections: Compounds, Expected Compounds, and Merged Features		
X Data	Specifies the principal component to plot on the <i>x</i> axis.		
	Selections: PC1, PC2, PC3, PC4, or PC5		
Y Data	Specifies the principal component to plot on the <i>y</i> axis.		
	Selections: PC1, PC2, PC3, PC4, or PC5		
Center and Scale	Centers and scales the data.		
Use Normalized Areas	Uses the normalized data.		
	Available if the processing workflow includes the Normalize Areas node.		
Scores Plot page	Displays a plot of one principal component versus a second principal component.		
Loadings Plot page	Displays the compounds in the selected table plotted against the selected principal components.		
	This plot is interactive with the selected results table.		
Variances Plot page	Displays the proportion and the cumulative proportion of the variance contributed by each principal component.		

Table 92. Principal Component Analysis parameters

Working with the Descriptive Statistics View

Use the Descriptive Statistics view to visually compare the statistics of the peak areas for all compounds currently displayed in the Compounds table or the Expected Compounds table as a box-and-whisker plot.

A box-and-whisker plot displays the data for a variable as a rectangular box with a set of whiskers at each end (Figure 103). The line through the rectangle represents the median value in the data set. The lower portion of the rectangle represents the data points that fall within the second quartile and the upper portion represents the data points that fall within the third quartile. The circles that fall outside the fence whiskers are outliers.

The application uses the following equations to calculate the upper and lower fences:

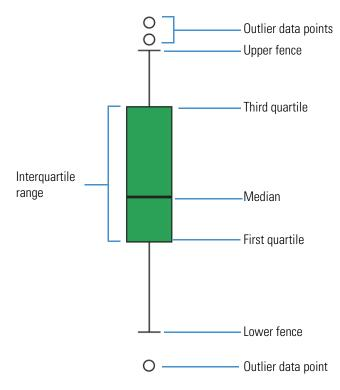
Interquartile range (IQR) = Quartile 3 (Q3) – Quartile 1 (Q1)

Upper fence = $Q3 + IQR \times 1.5$

Lower fence = $Q1 - IQR \times 1.5$

Note To calculate the quartiles, the application uses a method that is similar to the type 6 method in the R statistical computing software.

Figure 103. Box-and-whisker display of the data distribution



By default, the graph displays the data for all the samples and duplicates the grouping on the Grouping and Ratios page of the analysis. Each group appears in a different color. The legend shows the colors of the sample groups.

Use the Group By check boxes in the collapsible pane to change the grouping. Use the Filter By check boxes to remove samples from the plot.

For more information about the shortcut menu commands, see "Copying or Saving Graphical Views for Publication" on page 291.

Follow these procedures as needed:

- To copy or save the data as an image
- To copy or save the data as editable text

- To view the table row for an outlier data point
- To select the check box for an outlier data point
- To select the check boxes for all of the visible data points
- To export the outlier data points to a spreadsheet

To copy or save the data as an image

- 1. Right-click the plot and choose **Show Legend**.
- 2. Right-click the plot and choose Copy > Image to copy an image to the Clipboard.

-or-

Right-click the plot, choose **Export > Image As**, and select an image type to save the data to an image file.

* To copy or save the data as editable text

Right-click the plot and choose **Copy > Data** to copy the text to the Clipboard.

You can paste this text to Notepad, an Excel spreadsheet, and so on. The data appears in a columnar format.

-or-

Right-click the plot and choose **Export > Data As** to save the data to a text file.

The file contains two data sets. The first set consists of these columns from left to right: Groups, Name, Minimum Value, Maximum Value, Std. Deviation, Mean, Median, Q1 Value, Q2 Value, and Q3 Value. The second set lists the outlier data points and consists of these columns from left to right: Groups, Name, and Outlier.

✤ To view the table row for an outlier data point

In the plot, double-click the data point.

To select the check box for an outlier data point

In the plot, right-click the data point and choose Check Point.

In the result table, the check box is selected for the corresponding compound or expected compound.

* To select the check boxes for all of the visible data points

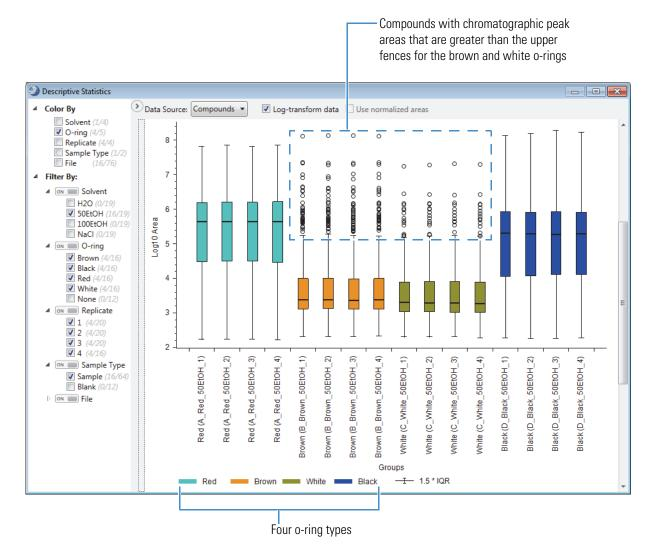
Right-click the plot and choose Check All Visible Points.

***** To export the outlier data points to a spreadsheet

1. In the view, zoom in on the outlier points so that they are the only visible points on the screen.

Figure 104 shows a box-and-whisker plot for the chromatographic peak areas of compounds that leach out of four different o-ring types. The compounds leached from the red and black o-rings show more variation; however, the peak areas for all of the these compounds fall within the lower and upper fences. The compounds leached from the brown and white o-rings show less variation (in peak area); however, the peak areas for numerous compounds fall outside the upper fences.

Figure 104. Box-and-whisker plot for the leachable compounds in four types of o-rings



- 2. Right-click the plot and choose **Check All Visible Points** to select the rows for these outliers in the Compounds table.
- 3. In the Compounds table, sort the checked rows by the variable of interest. For example, sort the rows by molecular weight, retention time, or both.
- 4. Right-click the Compounds table and choose **Export > Export to Excel**.
- 5. In the Export to Excel dialog box, select a folder and name the file, select the **Checked Items Only** check box, and click **Export**.

Table 93 describes the components of the Descriptive Statistics view.

Parameter	Description			
Data Source	Specifies the result table for the source data.			
	Selections: Compounds or Expected Compounds			
Log-transform Data	Determines whether the data appears in a linear scale or the \log_{10} scale.			
	Selecting this check box transforms the area counts to the \log_{10} scale.			
Use Normalized Areas	S Select to display normalized chromatographic peak areas.			
	Available when the processing workflow includes the Normalize Areas node.			
Graph				
x axis	Displays the name of the sample group.			
<i>y</i> axis	Displays the area in a linear scale or in a \log_{10} scale.			
Outlier data points	The circles represent outlier points. When the Show Position Tooltips command is enabled, placing the cursor over a data point displays the following information: MW, RT, Max. Area, and #Adducts.			

Table 93. Descriptive Statistics view parameters (Sheet 1 of 2)

Parameter	Description
Rectangle	The rectangles represent the second and third quartiles for the data set.
	When the Show Position Tooltips command is enabled, placing the cursor over the rectangle displays the following information: filename, group, maximum value (including the outliers), 3 rd quartile, median, 1 st quartile, minimum value (including the outliers).
Legend	By default, the legend is hidden.
	Choosing Show Legend from the shortcut menu displays the legend. The legend shows the group colors.

Table 93. Descriptive Statistics view parameters (Sheet 2 of 2)

Working with the Differential Analysis View

Use the Differential Analysis view to view a volcano plot of the differential analysis performed during data processing or to run a new differential analysis.

Follow these topics as needed:

- Reviewing the Initial Differential Analysis
- Changing the Analysis Settings for a Differential Analysis
- Running a New Differential Analysis
- Running a New Differential Analysis
- Differential Analysis View Parameters and Shortcut Menu Commands

Reviewing the Initial Differential Analysis

* To open the Differential Analysis view and review the initial analysis

- 1. Open the result file for an analysis that included at least two raw data files.
- 2. In the menu bar, choose **View > Differential Analysis**.

If the processing workflow included the Differential Analysis node, the Differential Analysis view opens with the differential analysis from data processing. The ratios in the Comparison list match the ratios on the Grouping and Ratios Summary page of the Summaries view. The initial p-value setting is $0.05 \ (-\log_{10} 0.05 = 1.3)$ and the initial \log_2 Fold change setting is 1 (a ratio of 2 to 1). Depending on the setting for the \log_{10} Transform parameter in the Differential Analysis node, the *y*-axis scale spans the p-value range (0-1) or the $-\log_{10}$ p-value range $(-\log_{10} 0 = \text{Infinity}, -\log_{10} 1 = 0)$.

3. To show the legend, right-click the plot and choose **Show Legend**.

The data points are color-coded, as defined by the legend.

Figure 105 shows a comparison between two extraction solvents—water and an aqueous solution of 50% ethanol.

Figure 105. Differential Analysis view with the analysis run during data processing



Changing the Analysis Settings for a Differential Analysis

To change the analysis displayed in the volcano plot

- Select a different ratio from the Comparison list. This changes the data points in the plot.
- Change the p-value setting by dragging the slider. This changes the *y*-axis range of the shaded areas.

Figure 106 shows the effect of changing the p-value setting.

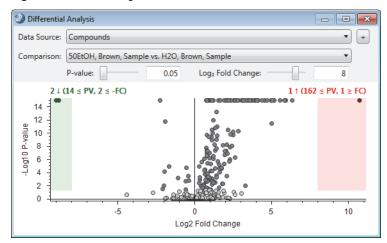
Differential Analysis - - -Data Source: Compounds • + Comparison: 50EtOH, Brown, Sample vs. H2O, Brown, Sample • Log₂ Fold Change: P-value: 1E-12 1 3 ↓ (3 ≤ PV, 19 ≤ -FC) 57 ± (66 < PV, 155 > FC) 14 12 -Log10 P-value 10 8 6 4 2 0 -5 5 10 0 Log2 Fold Change

Figure 106. P-value decreased from 0.05 to 1E–12

• Change the fold change setting by dragging the slider. This changes the *x*-axis range of the shaded areas.

Figure 107 shows the effect of changing the Log_2 Fold Change setting.

Figure 107. Fold change increased from 1 to 8



Running a New Differential Analysis

To run a new differential analysis

Note Use the Perform New ANOVA Calculation dialog box to set up and compare different ratios than those on the Grouping and Ratios page of the analysis.

1. In the Differential Analysis view, click the **Perform New ANOVA Calculation** icon, **±**], to the right of the Data Source box.

The Perform New ANOVA Calculation dialog box opens (Figure 108).

· · · · · · · · · · · · · · · · · · ·							
Differentia	l Analys	sis					
Data Source:	Comp	ounds					• +
Comparison:	50EtC	H, Brown, Samp	le vs. H2O, 0.05		iple Id Change:	Π	▼
9.1.1		e. V, 19 ≤ -FC)	0.05	LUYZTU	ia change.	114 + (1	.62 ≤ PV, 155 ≥ FC)
14		Perform Ne	w ANOVA	Calculation			•
12 - 10 - 10 - - - - - - - - - - - - - - - - - - -		Data Source: Fixed Effect:	Compound Solver	ds nt ate le Type		•	
0 1	· · ·	Reference:				•	10
0	N N S.g Sig		Transform o Use normal Use nested	ized areas			
•	Sig		Ca	lculate	Canc	el	

Figure 108. Perform New ANOVA Calculation dialog box

- 2. In the Data Source list, select a different result table if applicable.
- 3. In the Fixed Effect area, select one or more variables to create new sample groups.
- 4. In the Reference list, select one of the new sample groups as the denominator for the group ratios.
- 5. (Optional) To transform the data to the \log_{10} scale, select the **Transform Data** check box.
- 6. (Optional) To normalize the chromatographic peak areas, select the **Use Normalized Areas** check box. This check box is available if the processing workflow included the Normalize Areas node.
- 7. (Optional) To calculate p-values for a nested design, select the **Use Nested Design** check box. This check box is available if the original analysis included a biological study factor and nested sample groups.

8. Click Calculate.

The Perform New ANOVA Calculation dialog box closes and the Pending Results area appears at the bottom of the view. The Pending Results area displays the data source (result table), the reference group, and the progress of the new calculation.

Pending Results:		
Data Source	Reference	Progress
Compounds	100EtOH, Black	Cancel

When the calculation is complete, the new ratio (*sample group* vs. *reference group*) appears in the Comparison list.

9. To view the analysis, select the data source in the Data Source list and the new ratio in the Comparison list.

The volcano plot updates with the selected analysis.

Differential Analysis View Parameters and Shortcut Menu Commands

Table 94 describes the parameters in the Differential Analysis view.

Parameter	Description				
Data Source	Specifies the result table for the source data.				
	Selections: Compounds or Expected Compounds table				
Perform New ANOVA Calculation	Opens the Perform New ANOVA Calculation dialog box for setting up a new differential analysis.				
Comparison	Specifies the ratio for comparison.				
	Selections: Generated ratios on the Grouping and Ratios page of the analysis and any new analyses that you have run and saved				
	To save the differential analyses run in the Differential Analysis view, save the result file.				
P-value	Specifies the p-value for the null hypothesis.				
	Use the slider to change the value in the corresponding box.				
	Default: 0.05 (-log ₁₀ 0.05 = 1.3)				
Log ₂ Fold Change	Specifies the fold change (ratio in the log base 2 scale) between the sample group and the reference group. This value creates an upper and lower threshold for each group ratio.				
	Use the slider to change the value in the corresponding box. Data points that fall outside the upper and lower thresholds are in the shaded regions.				
	Default: 1 (two-fold change)				
Graph					
<i>x</i> axis	Displays the log ₂ fold change.				
<i>y</i> axis	Displays the p-value in a linear scale or in a –log ₁₀ scale.				
Analysis summaries	The summaries display the number of data points with a p-value above the statistical significance level and a fold change outside th empirical threshold.				

Table 94. Differential Analysis view parameters (Sheet 1 of 2)

Parameter	Description	
Shaded regions	The region shaded in red identifies the data points that are significantly different (populations differ based on the p-value setting) and that fall outside the lower fold change threshold.	
	The region shaded in green identifies the data points that are significantly different (populations differ based on the p-value setting) and that fall outside the upper fold change threshold.	
Legend	By default, the legend is hidden.	
	 Choosing Show Legend from the shortcut menu displays the legend. The legend contains color-coded circles for these conditions: Nonsignificant and does not meet FC (fold change) threshold Nonsignificant and greater than upper FC threshold Nonsignificant and less than lower FC threshold Significant and greater than upper FC threshold 	

Table 94. Differential Analysis view parameters (Sheet 2 of 2)

Table 95 describes the shortcut menu commands for the Differential Analysis view.

Command	Description			
Show Position Tooltips	Displays information about the compound.			
Zoom Out	Undoes the last zoom.			
Undo All Zoom/Pan	Zooms out to the full plot range.			
Copy and Export commands	See "Copying or Saving Graphical Views for Publication" on page 291.			
Show Legend	Displays the legend for the color-coded data points.			
source (the Compounds tab	points (circles) represent compounds in the selected data le or the Expected Compounds table). Selecting a data point check box for the corresponding compound in the data source			
Check Point	To activate this command, point to a data point.			
	Selects the data point in the view and the check box for the corresponding compound in the data source table.			
Uncheck Point	To activate this command, point to a selected compound.			
	Returns the data point to its original color clears the corresponding check box in the data source table.			

Table 95. Shortcut menu commands for the Differential Analysis view (Sheet 1 of 2)

Command	Description
Check All Visible Points	Selects all the visible data points.
Uncheck All Visible Points	Clears the check boxes for all the visible data points.
Check All Up-Regulated Points	Selects all the data points in the pink-shaded region of the plot.
Check All Down-Regulated Points	Selects all the data points in the green-shaded region of the plot.

Table 95. Shortcut menu commands for the Differential Analysis view (Sheet 2 of 2)

Working with the Partial Least Squares Discriminant Analysis View

Use the Partial Least Squares - Discriminant Analysis view to determine whether two groups are different and to identify the variables that contribute to that difference.

See these topics:

- Identifying a Set of Compounds to Discriminate Between Groups
- Partial Least Squares Discriminant Analysis View Parameters

Identifying a Set of Compounds to Discriminate Between Groups

- To determine a set of compounds that you can use to discriminate the experimental variables by
- 1. Open a result file from an analysis with study factors.
- 2. From the menu bar, choose View > Partial Least Squares Discriminant Analysis.

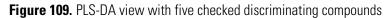
The Partial Least Squares Discriminant Analysis view opens to the right of the tabbed result tables.

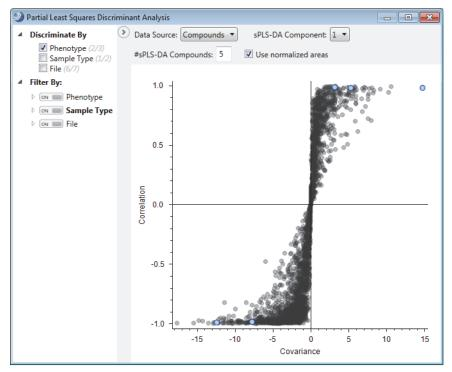
- 1. In the collapsible pane on the left, under Discriminate By, select the study factors that you want to discriminate by.
- 2. On the right above the plot, in the Data Source list, select the data source.
- 3. In the #sPLS-DA Compounds box, type the number of compounds that you want to use to differentiate the selected study factors.
- 4. To update the plot, click anywhere in the plot.

The orange circles represent the discriminating compounds.

5. Right-click the plot and choose Check All sPLS-DA Points.

The application selects the check boxes of the corresponding compounds in the selected data source. The blue circles represent the checked discriminating compounds (Figure 109).





6. To view the list of compounds that you can use to differentiate the selected experimental variables, set up a result filter to display only the checked compounds in the selected result table.

Partial Least Squares - Discriminant Analysis View Parameters

Table 96 describes the parameters in the Partial Least Squares - Discriminant Analysis view.

Parameter	Description		
Data Source	Specifies the result table for the source data.		
	Selections: Compounds or Expected Compounds table		
sPLS-DA Component	Specifies the analysis component.		
	Selection: 1 to 6		

Table 96. Partial Least Squares - Discriminant Analysis view parameters (Sheet 1 of 2)

Parameter	Description
#sPLS-DA Compounds	Specifies the number of compounds that when used together can discriminate between the values for the study variables selected under Discriminate By.
Use Normalized Areas	Uses the normalized areas in the result file.
	Available when the processing workflow includes the Normalize Areas node.

 Table 96.
 Partial Least Squares - Discriminant Analysis view parameters (Sheet 2 of 2)

Viewing KEGG Pathways

Use the KEGG Pathways view to view the reaction pathway of a detected compound in a KEGG pathway.

***** To view the KEGG pathways for a compound in the Compounds table

- 1. Open a result file that contains mapped KEGG pathways—that is, a result file from an analysis that included the Map to KEGG Pathways node.
- 2. In the main Compounds table, select a compound.
- 3. Below the main Compounds table, click **Show Related Tables** to show the related tables.
- 4. Click the **KEGG Pathways** tab to make it the active result table.
- 5. In the KEGG Pathways table, select the pathway that you want to view.
- 6. In the menu bar, choose View > KEGG Pathways.

The selected KEGG pathway opens. The blue circle indicates the selected compound in the KEGG pathway. The red circles indicate related compounds that were not found in the mzCloud database. The green circles indicate related compounds that were found in the mzCloud database. Figure 110 shows the selections of caffeine in the Compounds table and caffeine metabolism in the KEGG Pathways table. The related Compounds tables shows the compounds related to caffeine, which appear as red circles in the pathway diagram.

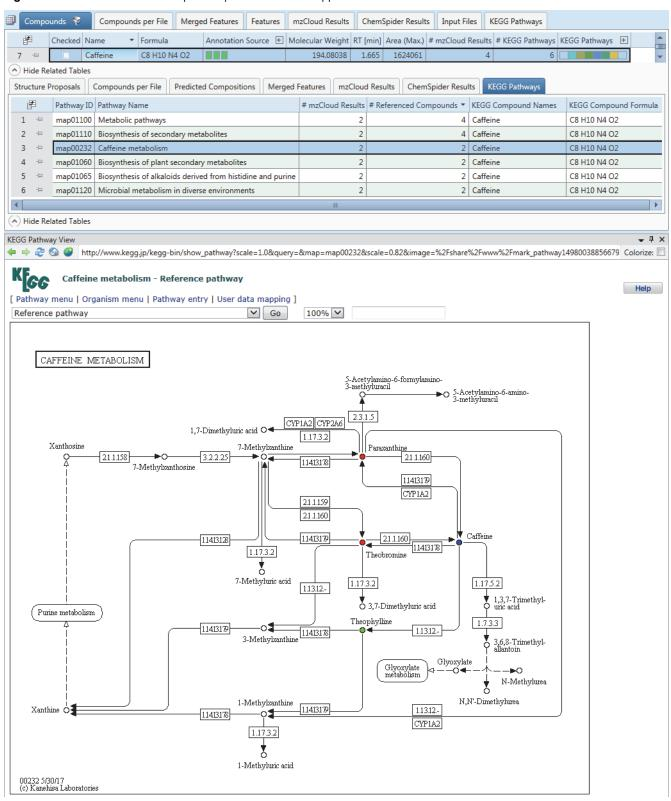


Figure 110. Caffeine metabolism pathway where caffeine appears as a blue circle

Viewing BioCyc Pathways

When the result file's processing workflow includes the Map to BioCyc Pathways Node, use the BioCyc Pathways view to view the mapped pathways for a selected compound.

* To view the BioCyc Pathways for a compound

- 1. Open a result file for an analysis that maps the detected compounds to their BioCyc pathways.
- 2. In the main table panel, click the **BioCyc Pathways** tab.
- 3. Select a pathway of interest.
- 4. Below the BioCyc Pathways table, click Show Related Tables.
- 5. Click the **BioCyc Results** tab.
- 6. In the BioCyc Results table, select a compound of interest (Figure 111).

BioCyc ×										
Comp	ounds	Compounds per File	e Featur	es BioCyc Results	Input Files	BioCyc Pathwa	ys			
Ē										
-	META:PWY-7300 ecdysone and 20-hydroxyecdysone bi									
2 +=	META:PWY-6061 bile acid biosynthesis, neutral pathway					10				
Hide R	Related Ta					- '				
Ē	Struct	ure		Name		Formula	Molecular Weight	BioCyc ID	BioCyc DB	
1 🗢	م	H H H H	ОН	3-dehydro-2,22-dideo	xyecdysone	C27 H42 O4	430.30831	CPD-15505	META	
2 눧	O;		Koh	3-dehydro-2-deoxyec	dysone	C27 H42 O5	446.30322	CPD-15506	META	
3 ⇔	HO	H OH H H OH H H	<u>ل</u>	2,22,25-trideoxyecdys	one	C27 H44 O3	416.32905	CPD-15499	META	
4 ⇔	но	H H H	Кон	2,22-dideoxyecdysone	2	C27 H44 O4	432.32396	CPD-15498	META	
5 ⇔	но	H H H H H H H H H H H H H H H H H H H	Ko#	3-dehydroecdysone		C27 H42 O6	462.29814	CPD-331	META	
6 ⇔	но	H H H H H H H H H H H H H H H H H H H	Ļан	2-deoxyecdysone		C27 H44 O5	448.31887	CPD-15497	META	
7 👳	0			3-dehydro-2,22,25-de	oxyecdysone	C27 H42 O3	414.31340	CPD-15500	META	

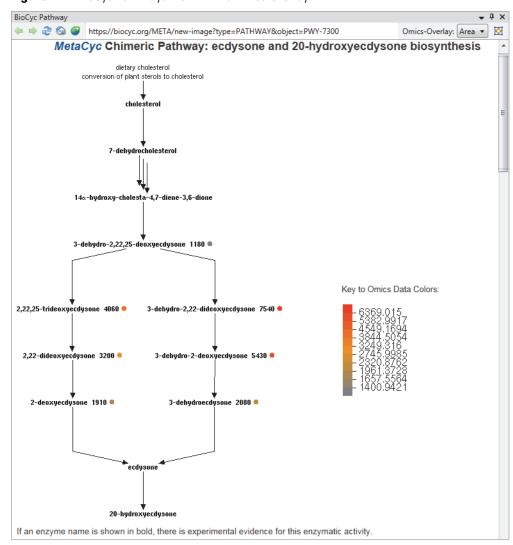
Figure 111. BioCyc Results table for the selected pathway

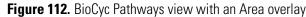
7. From the menu bar, choose **View > BioCyc Pathways**.

The BioCyc Pathways view opens to the right of the result tables.

- 8. In the upper right corner of the view, click the **Fit Pathway to Browser** icon, 🔯.
- 9. In the Omics-Overlay list (to the left of the icon), select Area.

Figure 112 shows a BioCyc pathway with an Area overlay.





Viewing Metabolika Pathways

Use the Metabolika Pathways view to view the mapped pathways for a selected compound.

To view a mapped pathway for a compound

- 1. Open a result file from an analysis that included the Map to Metabolika Pathways node.
- 2. In the Compounds table, select the compound.
- 3. Click Show Related Tables.
- 4. In the Related Tables panel, click the Metabolika Pathways tab.
- 5. Select the pathway that you want to view.

6. From the application menu bar, choose View > Metabolika Pathways.

The view opens to the right of the Compounds table and displays the selected pathway. The structure for the selected compound is blue, the structures for other detected compounds are red, and the structures for undetected compounds in the pathway are black.

7. To overlay a data source, select the source from the Overlay Data Source list and type or select a cell size from **6** to **30** pixels in width in the Overlay Cell Size box.

The data source selections depend on the processing workflow, but always include the annotation source, chromatographic peak area, or Metabolika pathways graphic.

Viewing the Corrected Retention Times of the Alignment Features

When the result file's processing workflow includes the Align Retention Times Node and the analysis includes more than one input file, use the Retention Time Correction view to inspect the regression curves for the alignment features.

To open the Retention Time Correction view

- 1. Open a result file for an analysis with multiple input files.
- 2. In the menu bar, choose **View > Retention Time Corrections**.

In the default layout, the Retention Time Correction view opens to the right of the tabbed result tables. Because you need to select a row in the Input Files table, the view is empty.

- 3. In the main result tables, click the Input Files tab.
- 4. Do the following as needed:
 - To view the regression curve for one input file as well as the prediction interval for the file's landmark features, select one input file (Figure 113).

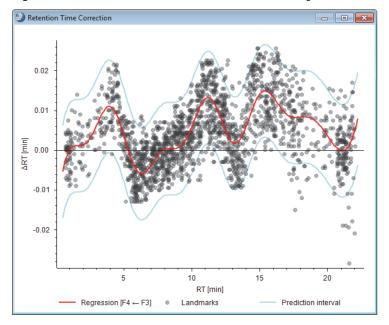


Figure 113. Retention Time Correction view with the regression curve for one input file

• To view overlaid regression curves for multiple input files, select multiple input files (Figure 114).

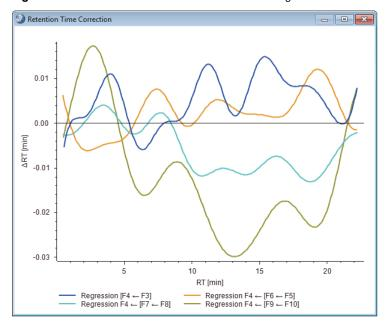


Figure 114. Retention Time Correction view with regression curves for multiple input files

Viewing the QC-Based Compound Area Correction Curves

When the result file's processing workflow includes the Normalize Areas Node and the analysis includes samples with the sample type assignment of Quality Control, use the Compound Area Corrections view to inspect the corrected peak areas for each detected compound in the Compounds Table across the input files.

For information about acquiring a set of raw data files with interspersed quality control samples, setting up the batch normalization criteria in the (post-acquisition) processing workflow, and setting up the Region and Language settings for the processing computer, see "Quality Control Samples for Batch Normalization" on page 34.

To open the Compound Area Correction view

- 1. Open a result file for an analysis with batch normalization—that is, and analysis that includes QC samples and a processing workflow with the Normalize Areas node.
- 2. In the Compounds table, select one or more compounds.
- 3. From the menu bar, choose View > Compound Area Corrections.

In the default layout, the Compound Area Correction view opens to the right of the tabbed result tables. The Compound Area Correction view displays a scatter plot of the areas for the selected compound on the y axis against the acquisition time for each input file on the x axis.

Figure 115 shows the Compound Area Corrections view for the compound selected in the Compounds table, with the following legend:

- Blue triangles represent the original compound areas for the QC samples.
- Orange triangles represent the corrected compound areas for the QC samples.
- Blue circles represent the original compound areas for the non-QC samples.
- Orange circles represent the corrected compound areas for the non-QC samples.
- The blue line is the regression curve for the original compound areas in the QC samples.
- The orange line is the regression curve for the corrected compound areas in the QC samples.

Pointing to a circle or triangle displays a tooltip with the input file name, the compound area, and the status—original or corrected.

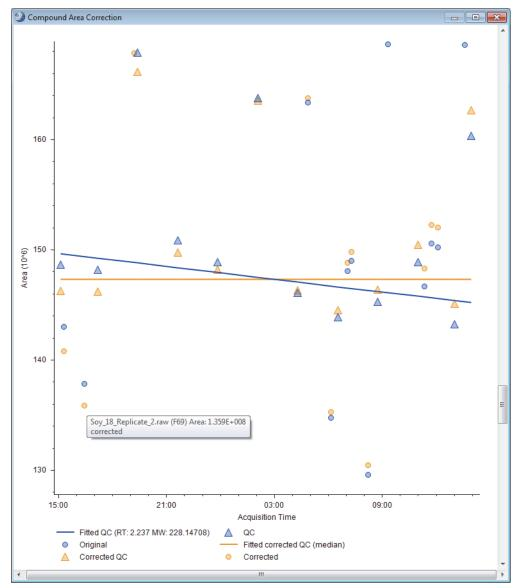


Figure 115. Compound Area Corrections view for a selected compound

Viewing the FISh Scoring Queue

Use the FISh Scoring Queue view to view the progress of FISh scoring.

- To open the FISh Scoring Queue view
- 1. Open a result file.
- 2. From the menu bar, choose **View > FISh Scoring Queue**.

The FISh Scoring Queue opens to the right of the result tables.

- 3. Start the FISh scoring process as follows:
 - a. Do one of the following for compounds with assigned structures:
 - In the Compounds table or the Expected Compounds table, right-click a compound and choose **Apply FISh Scoring**.
 - In any of the following related search results tables, right-click a compound and choose **Add to Structure Proposals and Apply FISh Scoring**.
 - In a Structure Proposals table, right-click a compound and choose either **Apply FISh Scoring to Selected** or **Apply FISh Scoring to All**.

The Settings dialog box opens.

b. In the Settings dialog box, make the appropriate selections and click OK.

-or-

- a. Do one of the following for compounds without assigned structures:
 - In the Compounds table or the Expected Compounds table, right-click a compound and choose **Edit Compound Annotations**.
 - In a Structure Proposals table, double-click a compound.

The Compound Annotation Editor dialog box opens.

- b. Enter the structure by opening a structure file, drawing the structure, or running a ChemSpider search.
- c. Click the FISh Scoring tab.
- d. Make the appropriate selections for the FISh scoring algorithm.
- e. At the bottom of the dialog box, select the **Apply FISh Scoring** check box and click **Save**.

Running a Hierarchical Clustering Analysis

Use the Hierarchical Cluster Analysis view to visualize the correlation between detected compounds and selected samples in a two-dimensional array of color-coded rectangles (heat map) where each rectangle represents the relative amount (by area) of a specific compound in a specific sample.

The application uses an agglomerative (bottom-up) approach to find the similarities between samples and compounds. Initially, the hierarchical cluster analysis assigns each compound to its own singleton cluster. The analysis then proceeds iteratively, at each stage joining the two most similar clusters into a new cluster, continuing until there is one overall cluster represented by a dendogram.

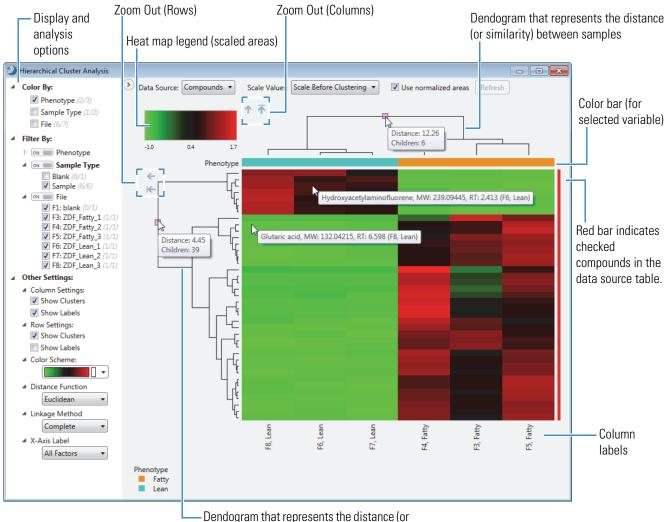
Figure 116 shows the results of a hierarchical cluster analysis for compounds detected in samples from six ZDF rats (phenotypes: Fatty and Lean).

The dendogram to the left of the heat map represents the distance (or dissimilarity) between the compound clusters. The width of each node is proportional to the distance (or dissimilarity) between two compounds at the lowest level or two compound clusters at the higher levels. The dendogram above the heat map represents the distance (or similarity) between samples. The height of each node is proportional to the distances (or similarity) between two samples at the lowest level or two sample clusters at the higher levels.

Pointing to a dendogram node displays the distance between clusters, and pointing to a heat map cell displays the compound and sample information.

By default, under Filter By, all the input files are selected, and under Other Settings, the Show Labels check box for the compound labels is clear.

Figure 116. Hierarchical cluster analysis for the compounds detected in six samples



dissimilarity) between compound clusters

To determine how to cluster the data, the application provides a choice of commonly used distance functions and linkage methods (see Table 98 and Table 99).

Note The Heat Map supports clustering for 2 to 2000 compounds.

To set up a hierarchical clustering analysis and work interactively with the heat map's cells, see these topics:

- Setting Up a Hierarchical Clustering Analysis
- Working Interactively with the Heat Map Grid
- Hierarchical Cluster Analysis View Parameters

Setting Up a Hierarchical Clustering Analysis

- * To set up the hierarchical clustering analysis
- 1. Filter the result table that you want to use as the data source. The analysis only uses the visible compounds in the table.

Note When you apply a filter that displays only the checked compounds in the selected result table, a red bar appears to the right of the heat map. If only a subset of the compounds is checked, the bar is discontinuous, with red indicating a checked compound and gray indicating an unchecked compound.

2. In the Data Source list, select the data source.

The Compounds table is the data source for untargeted analyses, and the Expected Compounds table is the data source for targeted analysis. For a combination analysis that generates a Compounds table and an Expected Compounds table, you can select either table as the data source.

- 3. In the Scale Values list, select whether to scale the heat map data before or after clustering or not at all.
- 4. To normalize the areas for the compounds, select the Use Normalized Areas check box.
- 5. In the left pane, under Filter By, select the check boxes for the samples to include in the analysis.

By default, all the sample files are selected.

Note By default, under Other Settings, the following check boxes are selected:

Column Settings: Show Clusters and Show Labels—Displays a dendogram above the heat map and sample labels below the heat map.

Row Settings: Show Clusters—Displays a dendogram to the left of the heat map.

- 6. In the left pane, under Other Settings, select these analysis options:
 - Under Distance Function, select how the analysis determines the distance between the data points, where the data points are files (columns) or compounds (rows) (Table 98).
 - Under Linkage Method, select how the analysis performs the clustering analysis (Table 99).
- 7. In the left pane, select these labeling options:
 - Under Color By, select whether to display color bars above the heat map to visually differentiate the samples by their study factor variables.
 - Under Other Settings > Row Settings, select whether to display the compound labels.

Selecting the Show Labels check box under Row Settings displays the compound labels to the right of the heat map.

- Under Other Settings > X-Axis Label, select the label for the sample columns.
- 8. To run the analysis, click **Refresh**.

Working Interactively with the Heat Map Grid

After you run a hierarchical heat map analysis, you can enlarge the heat map, zoom in and out on the heat map, and point to specific cells in the heat map for more information.

Task	Instructions
Enlarge the heat map.	To increase the heat map width, drag its left or right edge.
	To increase its height, drag its top or bottom edge.
Zoom in on an area of the heat map.	Drag the pointer across the rectangular area of the heat map.
Zoom out of an enlarged area of the heat map by using the shortcut menu commands.	Right-click the heat map and choose one of the following: Zoom Out Column, Zoom Out Row, Undo All Column Zoom, or Undo All Row Zoom.
Zoom out of an enlarged area of the heat map by using the zoom icons.	To zoom out of a row selection, click the Zoom Out Row icon, , or the Undo All Row Zoom icon, , below the heat map legend.
	To zoom out of a column selection, click the Zoom Out Column icon, 1 , or the Undo All Column Zoom icon, 7 , to the right of the heat map legend.

Task	Instructions
View information about a heat map cell.	Point to the cell. The tooltip displays the compound's name (if available) its MW, RT, file ID, and the study factor value (if available).
	Uric acid, MW: 168.02814, RT: 1.136 (F4, Fatty)
View the distance value for a	Point to the node.
dendogram node.	The tooltip displays the distance value.
	For an analysis where the chromatographic peak areas are scaled before clustering, the distance values are scaled.
	For an analysis where the chromatographic peaks areas are not scaled or are scaled after clustering, the distance values are not scaled.

Hierarchical Cluster Analysis View Parameters

Table 97 describes the parameters in the Hierarchical Clustering Analysis view.

Table 97. Hierarchica	al Clustering Analysis v	view parameters (Sheet 1 of 3)
-----------------------	--------------------------	--------------------------------

Parameter	Description
Button	
Refresh	Runs the hierarchical clustering analysis on the compounds displayed in the selected data source (result table) and the samples selected in the left pane under Filter By.
Parameters in the right pane	
Data Source	Specifies the result table for the source data.
	Selections: Compounds table or Expected Compounds table

Parameter	Description	
Scale Values	Specifies if and when to perform a z-score transformation on the data points:	
	• None—The application does not scale the data. The heat map legend displays a scale in area counts, and the dendogram nodes display the distance in area counts.	
	• Scale After Clustering— Applies a z-score transformation after performing the hierarchical clustering. The heat map legend displays the range of the scaled values, and the dendogram nodes display the distance in area counts.	
	• Scale Before Clustering—Applies a z-score transformation before performing the hierarchical clustering. The heat map legend displays the range of the scaled values, and the dendogram nodes display the scaled distance values.	
Use Normalized Areas	When selected, normalizes the chromatographic peak areas for the selected compounds before running the analysis.	
Parameters in the left pa	ne	
Color By		
Displays the selected co	lor bars for the selected variables above the heat map.	
Filter By		
•	c boxes—except for the Blank sample type—are selected. Clearing a corresponding item from the analysis.	
Other Settings		
Column Settings		
Show Clusters	When selected, displays the dendogram for the samples above the heat map; that is, this selection displays the dendogram for the items selected under Filter By.	
Show Labels	When selected, displays the labels across the bottom of the heat	
Show Labers	map for the items selected under Filter By. The X-Axis Label	
Row Settings	map for the items selected under Filter By. The X-Axis Label	

Table 97. Hierarchical Clustering Analysis view parameters (Sheet 2 of 3)

Parameter	Description
Color Schemes	Specifies the color scheme for the heat map.
	Default: Available values Example and Missing values []
	Selection:
	Available values Missing values
Distance Function	Specifies the distance function to use for calculating the distance
	between data points (see Table 98).
	Default: Euclidean
Linkage Method	Specifies the method to use for hierarchical clustering (see Table 99).
	Default: Complete
X-Axis Label	Specifies the labels for the sample columns.
	Selections: All Factors, Selected Factors, File ID, Full File Name, and Sample Name

Table 97. Hierarchical Clustering Analysis view parameters (Sheet 3 of 3)

Table 98. Distance functions (Sheet 1 of 2)

Distance function	Description
Euclidean	Computes the Euclidean distance between two data vectors, which is the geometric distance in the multidimensional space.
Manhattan	Computes the city-block (Manhattan) distance between two vectors. The Manhattan distance between two data items is the sum of the differences of their corresponding components. In most cases, the Manhattan distance measure yields results similar to the simple Euclidean distance. Note, however, that the effect of outliers is dampened, since they are not squared.
Maximum	Computes the maximum distance on any one of the dimensions between two vectors. Use this function to define two objects as different if they differ in any one of the dimensions.

Distance function	Description
Pearson	Computes the Pearson product-moment correlation, which is a measure for the shape similarity between two clusters.
Squared Euclidean	Computes the squared Euclidean distance between two data vectors. The Euclidean Squared distance metric uses the same equation as the Euclidean distance metric, but does not take the square root.

Table 98. Distance functions (Sheet 2 of 2)

Table 99. Linkage methods

Linkage method	Description
Average	Computes the distance between two clusters as the average distance between all pairs of objects in the two different clusters.
Centroid	Computes the distance between two clusters as the difference between centroids. The centroid of a cluster is the average point in the multidimensional space.
Complete	Computes the distance between two clusters as the greatest distance between any two objects in the different clusters (furthest neighbors).
Median	Computes the distance between two clusters as the difference between centroids, using the size of each cluster as a weighting factor.
Single	Computes the distance between two clusters as the distance of the two closest objects (nearest neighbors) in the clusters.
WARD	Computes the distance between two clusters using Ward's method, which uses an analysis of variance approach to evaluate the distances between clusters. The smaller the increase in the total within-group sum of squares as a result of joining two clusters, the closer they are. The within-group sum of squares of a cluster is defined as the sum of the squares of the distance between all objects in the cluster and the centroid of the cluster. Ward's method tends to produce compact groups of well-distributed size.
Weighted Average	Computes the distance between two clusters as the average distance between all pairs of objects in the two different clusters, using the size of each cluster as a weighting factor.

Running an mzLogic Analysis

You can use the mzLogic Analysis view in two ways:

- To run an mzLogic analysis for a data set that was processed with a workflow that did not include the Apply mzLogic node
- To update an older analysis with new similarity results from the mzCloud spectral database

Note The application does not save the results from this mzLogic analysis to the result file; however, you can add suitable candidates to the Structure Proposals table for a compound and apply FISh scoring.

The ranking score provided by an mzLogic analysis is not a probability score. It is only a measure of how similar a putative structure is to closely matching structures in the mzCloud spectral database.

To run and review an mzLogic Analysis, see these topics:

- Performing an mzLogic Analysis
- Reviewing the Results of an mzLogic Analysis
- mzLogic Analysis View Parameters

Performing an mzLogic Analysis

To perform an mzLogic analysis

- 1. Open a result file from an untargeted analysis.
- 2. Select a compound in the Compounds table.
- 3. From the application menu bar, choose **View > mzLogic Analysis**.

The mzLogic Analysis view opens to the right of the Compounds table. If the selected result file does not include results from an identity search, the Candidates area is empty. Otherwise, the Candidates area contains candidates from the identity search nodes.

- 4. To add candidates from the ChemSpider database, click ChemSpider and run a search.
- 5. To run forward and reverse similarity searches on the candidates, click **mzCloud Similarity**.

If the application finds similar structures in the mzCloud database, it populates the Similar Structures from mzCloud area and the Calculate Score button becomes available. The structure tiles in this area include an mzCloud Match score and a Coverage value at the top and a formula and delta mass value at the bottom.

6. To display an mzLogic score for each candidate, click Calculate Score.

Reviewing the Results of an mzLogic Analysis

* To review the results of an mzLogic analysis

- 1. In the mzLogic Analysis view, open the Similar Structures from mzCloud area.
- 2. In the Candidates area, click a candidate to select it.

In the Similar Structures form mzCloud area, the matching portions of the similar structures are highlighted in blue.

mzLogic Analysis View Parameters

Parameter	Description
ChemSpider	Opens the ChemSpider Search dialog box (see "Finding a Structure in the ChemSpider Database" on page 541).
mzCloud Similarity	Runs a forward and reverse mzCloud similarity search.
Calculate Score	Calculates the mzLogic scores for the available candidates.
	Available when the Candidates and Similar Structures from mzCloud areas are populated.
Back	Returns the focus to the previously selected compound in the Compounds table.
	Available when the Auto Refresh check box is clear and you select a different compound in the Compounds table.
Refresh	Refreshes the Candidates area when you select another compound.
	Available when the Auto refresh check box is clear.
Auto refresh	Select so that when you select another compound, the application automatically refreshes the Candidates area.
Size	Controls the size of the structure tiles in the Candidates area.
	Default: Medium
	Selection: Small, Medium, or Large

Table 100.mzLogic Analysis view parameters (Sheet 1 of 2)

Parameter	Description
Candidates area	Displays structure candidates as tiles. The structure candidates are provided by identity searches during data processing or a separate ChemSpider search from the result file. If the identity searches find duplicate structures, the application consolidates them.
	Right-clicking a tile displays a shortcut menu with the following commands:
	• Add to Structure Proposals—Adds the selected structure to the Structure Proposals table for the selected compound.
	• Add to Structure Proposals and Apply FISh Scoring—Opens the Settings dialog box for applying the FISh scoring algorithm. Then, runs the FISh scoring algorithm and adds the structure and FISh Coverage score to the Structure Proposals table for the selected compound.
Similar Structures from mzCloud area	Displays the results of the mzCloud similarity search for the selected compound.
	Selecting a candidate highlights the maximum common substructure in blue for the similar structures.
Sort By	Sorts the similar structures by the selected parameter.
	Default: mzCloud Match
	Selection: mzCloud Match, Forward Coverage, or Reverse Coverage

Table 100.mzLogic Analysis view parameters (Sheet 2 of 2)

Working with the Isotopologues Distribution Chart

If the analysis included the Analyze Labeled Compounds node, labeled samples, and at least one unlabeled reference sample, you can view a distribution chart of the detected isotopologues.

* To open the Isotopologues Distribution Chart and review the isotopologues

- 1. Open a result file for a stable isotope labeling analysis.
- 2. From the application menu bar, do one of the following:
 - Choose Window > Apply Layout > Stable Isotope Labeling.

The Isotopologues Distribution Chart, Trend Chart, and Metabolika Pathways view open as a tabbed group to the right of the result tables. Under Group By, the File check box is selected. -or-

• Choose View > Isotopologues Distribution Chart.

The Isotopologues Distribution Chart opens to the right of the result tables. Under Group By, all the check boxes are clear.

The compound's name, elemental composition, MW, and RT appear above the graph (Figure 117). The bars are colorized by input file.

Figure 117. Isotopologues Distribution Chart with no grouping selected



- 3. Do any of the following:
 - To display a tooltip for a bar, point to the bar.

Tip By default, the Show Position Tooltips feature is turned on. If a tooltip does not appear, click the chart. If a tooltip still does not appear, right-click the chart and choose **Show Position Tooltips**.

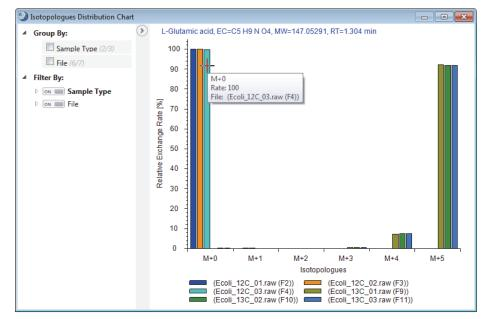
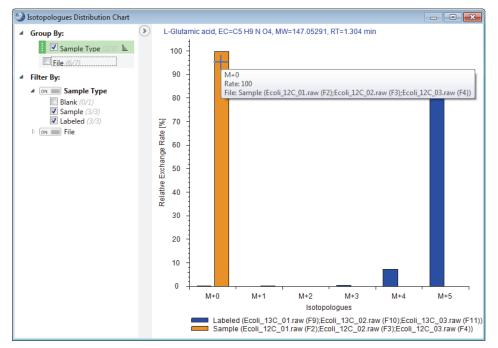


Figure 118. Isotopologues Distribution Chart with a tooltip

- To group the isotopologues by input file, under Group By, select the File check box.
- To group the isotopologues by sample type, under Group By, select the **Sample Type** check box (Figure 119).

Figure 119. Isotopologues grouped by sample type



9 Working with the Graphical Result Views Working with the Isotopologues Distribution Chart



Reviewing the Tabular Result Data

For information about the tabular data in the result files, see these topics.

Contents

- Common Result Tables
- Expected Compounds Result Tables
- Compound Detection Result Tables
- Compound Identification Result Tables
- Pathway Mapping Result Tables
- Compound Scoring Tables
- Differential Analysis Columns
- Descriptive Statistics Columns

For general information about opening result files and working with the result tables, see Chapter 8, "Reviewing the Analysis Results."

Common Result Tables

You can run a targeted analysis that finds expected compounds, an untargeted analysis that detects unknown compounds, or both types of analyses with the Compound Discoverer application.

For information about the result tables that are independent of the analysis type, see these topics:

- Adducts Table
- Chromatogram Peaks Table
- File Alignments Table
- FISh Trace Fragments Table

- Input Files Table
- Manual Peaks Table
- Merged Features Table
- Specialized Traces Table
- Structure Proposals Table

Adducts Table

Use the Adducts table to view the list of adducts in the Adducts library. By default, the Adducts table is hidden (see "Showing or Hiding Result Tables" on page 257).

Chromatogram Peaks Table

Use the Chromatogram Peaks table to view information about the quality of the chromatographic peak for the detected or expected feature. The Chromatogram Peaks table is related to the Features and Expected Features tables.

Table 101 describes the columns in the Chromatogram Peaks table.

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Area	Displays the integrated peak area.
Apex m/z	Displays the <i>m/z</i> value of the mass spectral peak at the chromatogram peak apex.
Apex RT [min]	Displays the retention time of the chromatogram peak apex.
Apex Intensity	Displays the intensity at the chromatogram peak apex.
Left RT [min]	Displays the start point for the chromatographic peak.
Right RT [min]	Displays the end point for the chromatographic peak.

 Table 101.Chromatogram Peaks table (Sheet 1 of 2)

Column	Description
Isotope Number (hidden)	(For chromatographic peaks detected by the Detect Compounds node) Displays the index number for the isotopic mass spectrum peak that the application used to create the XIC. The Detect Compounds node creates an XIC trace for each isotope.
	(For chromatographic peaks found by the Find Expected Compounds node) Always displays a value of 0, as the Find Expected Compounds node creates only one filtered XIC trace for each ion. The Find Expected Compounds node creates the filtered XIC trace by summing the intensity of all the mass spectrum peaks that match the theoretical isotope pattern. If even one required isotope is missing, the intensity of the XIC drops down to 0.
Peak Model	Displays the peak model for the chromatographic peak and includes information about the peak's width and symmetry.

Table 101.Chromatogram Peaks table (Sheet 2 of 2)

File Alignments Table

Use the File Alignments table to check the alignment process for each input file.

The Align Retention Times Node creates the File Alignments table.

Table 102 describes the columns in the File Alignments table.

Table 102	File Alignments table
-----------	-----------------------

Column	Description
Study File ID	Displays the study file ID of the sample where the algorithm has corrected the measured retention time of the detected features against a set of features from the reference file.
Ref. File ID	Displays the study file ID of the sample being used as a reference.
Kind	Displays the parameter that the algorithm used in the regression model. Retention time (RT) is the current regression model.
Description	Displays a description of the alignment process.
#Landmarks	Displays the number of features that the analysis used to align the specified file to the reference file.
Alignment Details	Displays additional information about the alignment.
RMSE	Displays the estimated error for the corrected retention times of the features in the selected input file as the root mean square error.

FISh Trace Fragments Table

Use the FISh Trace Fragments table to view the structures and the summed intensities of the expected fragment ions.

The Create FISh Trace Node creates the FISh Trace Fragments table when the Individual Traces parameter for this node is set to True.

Table 103 describes the columns in the FISh Trace Fragments table.

Table 103.FISh Trace Fragments table

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Parent Compound	Displays the selected compound in the Create FISh Traces node.
Formula	Displays the elemental composition of the fragment ion.
Ion	Displays the ion description.
m/z	Displays the ion's mass-to-charge ratio.
TIC	Displays the total ion current for the fragment ion.
Mode	Displays the fragmentation mode.
File ID	Displays the integer that the application assigned to the input file.
Study File ID	Displays the study file ID of the input file.
Structure	Displays the ion's molecular structure.

Input Files Table

Use the Input Files table, which is common to all processing workflows, to view information about the input file set (Xcalibur RAW files) that the application processed to produce the result file. For a result file from an analysis of multiple input files, use the related File Alignments Table to check the chromatographic alignment of the features in the input files.

* To check the chromatographic alignment of the input files

- 1. Click the Input Files tab.
- 2. Select a row (input file).
- 3. Click Hide Related Tables below the main tables.
- 4. Click the File Alignments tab.
- 5. From the menu bar, choose View > Retention Time Corrections.

The Retention Time Corrections view displays the regression curve for the selected input file.

Table 104 describes the information displayed in the Input Files table.

Column	Description
Study File ID	Displays the identification number assigned by the Compound Discoverer application.
File Name	Displays the file name of the input file.
Creation Date	Displays the acquisition time stamp from the data system.
RT Range [min]	Displays the data acquisition time for the raw data file.
Instrument Name	Displays the mass spectrometer type used to acquire the raw data file.
Software Revision	Displays the software version of the instrument control software used to acquire the raw data file.
Instrument Hardware (hidden)	Displays the hardware version of the Thermo Scientific mass spectrometer or analog detector used to acquire the raw data file.
Ref. File ID	Displays the reference file that the retention time alignment algorithm used to align the landmark features.
Sample Type	Displays the sample type.
Study Factors	Displays the study factor value (item).

Manual Peaks Table

Use the Manual Peaks table to view information about the manual peaks that you add to the result file.

The Manual Peaks table is a main table. It is also related to the main Specialized Traces table. To create this table, you must add a manual peak to a specialized trace (see "To manually integrate chromatographic peaks" on page 299).

Table 105 describes the columns in the Manual Peaks table.

 Table 105.
 Manual Peaks table

Column	Description
Trace Type	Displays the trace type. The trace type can be any of the specialized traces, including Analog, UV, PDA, TIC, BPC, XIC, Pattern Trace, or FISh Trace.
Area	Displays the chromatographic peak area.
Left RT [min]	Displays the start point of the chromatographic peak.
Right RT [min]	Displays the end point of the chromatographic peak.
Study File ID	Displays the study file ID of the input file (see "Removing Input Files or Updating Their Location" on page 105).

Merged Features Table

Use the Merged Features table to view ion conflicts between the Features and Expected Features tables. Also use the Merged Features table to correlate the chromatographic peaks detected by an analog detector to the chromatographic peaks detected by the mass spectrometer.

The Merge Features Node adds the Merged Features table to the result file. The Merged Features table has the following primary related tables: Expected Compounds, Expected Compounds per File, Compounds, Compounds per File, and Manual Peaks. Figure 7 on page 23 shows the hierarchy of the result tables for a targeted analysis, and Figure 9 on page 27 shows the hierarchy of the result tables for an untargeted analysis.

Tip When you add the Merge Features node to the processing workflow, the application automatically connects the Find Expected Compounds node and the Detect Compounds node to it. The Merge Features node consolidates the chromatographic peaks detected by these two nodes.

Table 106 describes the columns in the Merged Features table.

Table 106. Merged Features table (Sheet 1 of 2)

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Name	Use this column to describe or name the found chromatographic peak. To name the chromatographic peak, type alphanumeric text in the Name table cell.
Apex m/z	Displays the area weighted average mass of all related features.
	$\frac{\sum m/z \times \text{Area}}{\sum m/z}$
RT [min]	Displays the area weighted average retention time of all related features (same $m/z \times RT$ dimensions within the specified tolerances).
	$\mathbf{\nabla}$ RT × Area
	$\frac{\sum RT \times Area}{\sum RT}$
	$\sum RT$
Max. Area	Displays the area of the largest chromatographic peak found in the data set for the current $m/z \times RT$ dimensions.
Ion Conflict	Indicates whether there is a conflict between the Detect Compounds and Find Expected Compounds nodes.
(■) Green	No conflict—Both nodes assigned the same ion to this chromatographic peak.
(]) Orange	Missing unknown component—Not found by the Detect Compounds node.
	Indicates that the expected compound might be in doubt because the Detect Compounds node did not detect it.
(□) Gray	No data to compare—The related node did not detect this ion.
(]) Red	Either the two nodes assigned different ions or one of the nodes assigned more than one ion to this m/z value and retention time.
Detect Compounds	Indicates whether the Detect Compounds node found the current feature in each input file.
(□) Gray	No matches found
(D) Green	Single match found
(D) Red	Multiple matches found

Column	Description
Find Expected Compounds	Indicates whether the Find Expected Compounds node found the current feature in each input file.
(□) Gray	No matches found
(D) Green	Single match found
(]) Red	Multiple matches found
Max. Areas (for each input file)	For each input file, displays the maximum chromatographic peak area for the features with the same $m/z \times RT$ dimensions (within the specified tolerances) found by the Find Expected Compounds node, the Detect Compounds node, or both nodes.
Note The Differential Analysis node generates these columns: Group Areas, Ratio, and Log2 Fold Change. For information about these columns, see "Differential Analysis Columns" on page 432.	
Comments	Use this column to store comments about the current feature (unique $m/z \times RT$ dimensions). This column accepts alphanumeric text and special characters.

 Table 106.Merged Features table (Sheet 2 of 2)

Specialized Traces Table

Use the Specialized Traces table to view traces created by the tracer nodes.

The Specialized Traces table lists the specialized traces that you requested in the processing workflow. For information about manually integrating chromatographic peaks in a specialized trace, see "To manually integrate chromatographic peaks" on page 299.

For information about the Chromatograms view, see "Working with the Chromatograms View" on page 294.

✤ To view a trace in the Chromatograms view

Select the trace of interest in the Specialized Traces table.

Table 107 describes the columns in the Specialized Traces table.

 Table 107.Specialized Traces table (Sheet 1 of 2)

Column	Description
Checked	Use this column to select the rows that you want to display in the results table and in reports after you apply result filters.
Study File ID	Displays the study file ID of the input file (see "Removing Input Files or Updating Their Location" on page 105.

Column	Description
Trace/Detector Type	 Displays the trace type generated during data processing. Create Mass Trace Node—TIC, BPC, or XIC Create Analog Trace Node—UV Trace, Total Scan, Spectrum Maximum, <i>Wavelength–Wavelength</i>, or Analog trace Create FISh Trace Node—FISh Trace Create Pattern Trace Node—Pattern Trace
Custom Label	Displays the text that you entered in the Custom Label box for the processing workflow node that generated the trace. You can edit the text in this column.
Description	Displays a description of the trace.
Spectrum File	Displays the file name of the raw data file that includes the trace.

 Table 107.Specialized Traces table (Sheet 2 of 2)

Structure Proposals Table

Use the Structure Proposals table to store custom structure proposals for the selected compound or expected compound in the result file.

For information about adding structure proposals to the table, see "Adding and Deleting Proposed Structures for a Compound" on page 261.

Table 108 describes the columns in the Structure Proposals table.

 Table 108.Structure Proposals table (Sheet 1 of 2)

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Structure	Displays the structure of the compound (see "Modifying the Result Page Layout" on page 253).
Name	Displays an application-generated name or a user-specified name for the compound.
	To edit this entry, click the table cell and type text in the box.
Formula	Displays the elemental composition formula of the compound.
Molecular Weight	Displays the molecular weight of the compound.

Column	Description
FISh Coverage	Displays the FISh coverage score that is based on the proposed structure (see "FISh Scoring for Targeted Compounds and Proposed Structures" on page 36).
Comments	Displays an application-generated comment or a user-specified comment for the compound.
	To edit this entry, click the table cell and type text in the box.

Table 108.Structure Proposals table (Sheet 2 of 2)

Expected Compounds Result Tables

For information about the result tables for a targeted workflow with the Find Expected Compounds node, see these topics:

- Expected Compounds Table
- Expected Compounds per File Table
- Expected Features Table
- Expected Formulas Table
- Related Structures Table
- Transformations Table

Expected Compounds Table

Use the Expected Compounds table to view information about the targeted compounds that the analysis finds. This result table contains all of the expected compounds that the analysis found in the input file set. The compounds are grouped by formula, molecular weight, and retention time.

The Group Expected Compounds Node adds the Expected Compounds table to the result file. The Descriptive Statistics post-processing node adds the descriptive statistics columns, and the Differential Analysis post-processing node adds the differential analysis columns (Ratio, Log2 Fold Change, P-value, and Adjusted P-value).

Table 109 describes the columns in the Expected Compounds table.

Table 109. Expected Compounds table (Sheet 1 of 3)

Column	Description
Checked	Use this column to select the rows that you want to display in the results table and in reports after you apply result filters.
Structure (hidden)	Displays the structure of the compound if available.

Column	Description
Name	Displays the user-specified compound name.
	To populate this cell, you can type a name or use the Edit Compound Annotation command.
Parent Compound	Displays the selected compound for each Generate Expected Compounds node that you connected to the Find Expected Compounds node. Before you can select a specific compound in the Generate Expected Compounds node, you must first add it to the compound library (see "Importing, Exporting, and Deleting Compounds" on page 498).
Formula	Displays the elemental formula of the parent compound and the theoretical formulas for the dealkylation and transformation products.
Molecular Weight	Displays the molecular weight (MW) of the compound calculated from the formula (neutral mass). Expected compounds include the parent compounds and their theoretical dealkylation and transformation products.
Dealkylated	When this column contains an X, the expected compound is the product of a dealkylation reaction.
Transformations	Displays the chemical transformation for the expected compounds that have undergone any of the user-specified transformations.
Composition Change	Displays the composition change caused by any dealkylation or dearylation reaction, any of the user-specified transformation reactions, or both.
RT (min)	Displays the weighted average of the retention times for this chromatographic peak in the input file set.
RT Tolerance (min.)	Displays the retention time tolerance setting in the Group Expected Compounds Node.
FISh Coverage	Displays the FISh coverage score from the FISh Scoring Node (see "FISh Scoring for Targeted Compounds and Proposed Structures" on page 36).
Area (Max.)	Displays the area of the largest chromatographic peak found for this compound (Formula × MW× RT) in the sample set.
#Adducts (hidden)	Displays the number of adducts found for this compound.
#MS2 (hidden)	Displays the number of available MS2 scans for the compound. In the MS2 scans are not from an Identification Only sample, you can view them by selecting the rows in the related Expected Features table.

Table 109. Expected Compounds table (Sheet 2 of 3)

Column	Description
Area (per input file)	Displays the area for the compound (Formula × MW× RT) in each sample (input file).
	To display the sample names, click the expand icon to the right of the column heading.
Note See the following the hidden descriptive	ng topics for information about the differential analysis columns and e statistics columns.
• Saa "Descriptive	Statistics Columns" on page 425 for information about the following

Table 109. Expected Compounds table (Sheet 3 of 3)

- See "Descriptive Statistics Columns" on page 435 for information about the following columns: Min. Area, Median Area, Mean Area, Q1 Area, Q3 Area, Area CV%, and Area SD.
- See "Differential Analysis Columns" on page 432 for information about the following columns: Group Areas, Ratio, and Log2 Fold Change.

Expected Compounds per File Table

Use the Expected Compounds per File table to review the expected compounds found in each input file. The uniqueness for each row is defined by the following expression:

Parent Compound × Formula × MW × Dealkylations and Transformations × RT

Clicking a row in the Expected Compounds per File table displays an XIC trace for the selected compound (Figure 120). The XIC trace is a summation of the related ion traces. The integrated peak area is shaded, the vertical red line indicates the chromatographic peak apex, and the triangle indicates the data point that corresponds to the retention time (RT) label. If the data has been chromatographically aligned (by using the Align Retention Times node), the RT values for the labeled data points might differ slightly from those for the closest MS1 scan in the spectral tree.

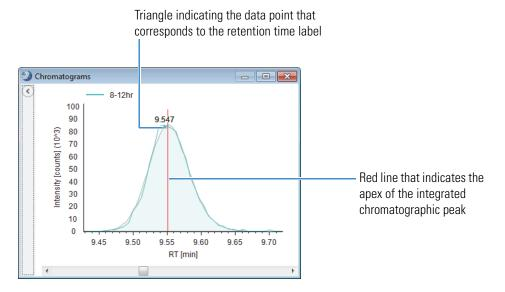


Figure 120. Expected compound trace for a single input file

The Find Expected Compounds Node creates the Expected Compounds per File table. The primary tables related to this table are as follows: Expected Compounds, Input Files, Expected Formulas, Merged Features, Expected Features, and Related Structures (see Figure 7 on page 23).

Table 110 describes the columns in the Expected Compound per File table. By default, some of these columns are hidden.

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Parent Compound	Displays the user-specified parent compound or compounds.
	You specify the parent compound or compounds in the Generate Expected Compounds node or nodes.
Formula	Displays the elemental formula of the expected compound.
Molecular Weight	Displays the molecular weight (MW) of the expected compound.
Dealkylated	Displays an X if the parent compound has undergone a dealkylation reaction.
Transformations	Displays the chemical transformations for the expected compound.
Composition Change	Displays the composition change caused by a dealkylation reaction, any of the user-specified transformation reactions, or both.

Table 110.Expected Compounds per File table (Sheet 1 of 3)

Column	Description
RT [min]	Displays the apex retention time (in minutes) of the largest chromatographic peak that the node found for the expected compound.
	The chromatographic peak is a composite peak of all of the ionic species (adduct ions) that the analysis found for the expected compound. The chromatographic peak area is the summed area of the adduct peaks.
FWHM [min]	Displays the width of the chromatographic peak at its half-height. Use this value to determine the best RT tolerance for peak grouping.
Best SFit [%]	Displays the best spectral fit value for the set of expected compound ions for the expected compound. The spectral fit value increases as the number of matching isotopes increases.
Best SD (hidden)	Displays the best spectral distance value between the theoretical and measured isotope pattern.
Max. #MI	Displays the maximum number of matching isotopes for any of the expected compound ions.
#Adducts	Displays the number of detected adducts. The analysis only detects the adduct ions that you specified for the Ions parameter in the Generate Expected Compounds node. With the default setting of [M+H]+1 only, the application finds only this one adduct ion species for each compound.
Area	Displays the summed chromatographic peak area for all of the expected compound ions (adducts) that make up the chromatographic peak.
	To display the table of expected compound ions for the expected compound, show the related tables and click the Expected Features tab.
Parent Area [%]	Displays the area of the selected component (Parent Compound × Formula × MW × Dealkylations and Transformations × RT) as a percentage of the total chromatographic peak area for the related components (Parent Compound × Formula × MW × Dealkylations and Transformations).

Table 110. Expected Compounds per File table (Sheet 2 of 3)

Column	Description
File ID	Displays the integer that the application assigned to the input file.
Study File ID	Displays the study file ID of the input file (see "Removing Input Files or Updating Their Location" on page 105).
	You can filter the data by using this integer; for example, the following filter reduces the table to the expected compounds found in one input file: Study File ID is equal to F1, F2, F3F11, and so on.

Table 110.Expected Compounds per File table (Sheet 3 of 3)

Expected Features Table

Use the Expected Features table to review the expected features (chromatographic peaks with the same $m/z \times RT$ dimensions) found across the input file set.

The Find Expected Compounds node creates the Expected Features table.

Table 111 describes the columns in the Expected Features table.

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Ion	Displays the ionized form of the compound.
Charge	Displays the charge of the ion.
Molecular Weight	Displays the molecular weight of the monoisotopic neutral compound.
m/z	Displays the mass-to-charge ratio of the ion.
ΔMass [Da]	Displays the mass difference, in daltons, between the theoretical mass of the ion and the measured mass.
ΔMass [ppm]	Displays the mass difference, in ppm, between the theoretical mass of the ion and the measured mass.
RT [min]	Displays the chromatographic retention time of the ion.
FWHM [min]	Displays the width of the chromatographic peak at its half-height in minutes.
SFit [%]	Displays the similarity score between the theoretical and measured isotope patterns as a percentage.
SD	Displays the spectral distance score.
#MI	Displays the number of matched isotopes for the ion.

Table 111. Expected Features table (Sheet 1 of 2)

Column	Description
Intensity	Displays the maximum intensity of all the related peaks per input file.
Area	Displays the summed area of all the related peaks (same expected compound) in the current input file.
Parent Area [%]	Displays the chromatographic peak area of the current ion as a percentage of the total chromatographic peak area at the current retention time (within the specified tolerance) for the expected compound selected in the Expected Compounds per File table.
File ID	Displays the integer that the application assigned to the input file.
Study File ID	Displays the study file ID of the input file (see "Removing Input Files or Updating Their Location" on page 105).

Table 111.Expected Features table (Sheet 2 of 2)

Expected Formulas Table

Use the Expected Formulas table to review the chemical formulas that the analysis found across the input file set. This table lists the theoretical compounds that the Generate Expected Compounds nodes predict by evaluating the effect of the user-specified dealkylation, dearylation, and transformation reactions on the user-specified parent compounds.

Clicking a row in the Expected Formulas table displays overlaid XIC traces for the selected expected compound, with one XIC trace for each input file where the compound is detected. Each XIC trace is a summation of the ion traces for the same neutral elemental composition (same molecular weight). By default, the Chromatograms view zooms in on the *x*-axis range of the detected peaks for the same expected compound.

The Find Expected Compounds Node creates the Expected Formulas table, which has the following primary related tables: Expected Compounds, Input Files, Expected Compounds per File, Related Structures, and Transformations.

For information about targeted processing workflows, see "Targeted Workflows for Expected Compounds" on page 16.

Table 112 describes the columns in the Expected Formulas table.

Column	Description
Checked	Use this column to select the rows that you want to display in the results table and in reports after you apply result filters.
Order (related table)	When you select the Expected Formulas table that is related to the Transformations table, this column displays the order of the selected transformation.

Table 112. Expected Formulas table (Sheet 1 of 2)

Column	Description
Parent Compound	Displays the selected compound for each Generate Expected Compounds node that you connected to the Find Expected Compounds node. Before you can select a specific compound in the Generate Expected Compounds node, you must first add it to the compound library (see "Importing, Exporting, and Deleting Compounds" on page 498).
Formula	Displays the elemental formula of the parent compounds and the theoretical formulas for the dealkylation and transformation products.
Molecular Weight	Displays the molecular weight (MW) of the expected compound. Expected compounds include the parent compounds and their theoretical dealkylation and transformation products.
Dealkylated	When this column contains an X, the expected compound is the product of a dealkylation reaction.
Transformations	Displays the chemical transformation for the expected compounds that have undergone any of the user-specified transformations.
Composition Change	Displays the composition change caused by any dealkylation or dearylation reaction, any of the user-specified transformation reactions, or both.
Area (Max.)	Displays the maximum summed chromatographic peak area for the expected formula in one of the input files.
	• When the result file contains data from only one input file, this area matches the summed chromatographic peak area for the expected formula.
	• When the result file contains data from more than one input file, this area comes from the input file with the largest summed chromatographic peak area for the expected formula

Table 112. Expected Formulas table (Sheet 2 of 2)

Related Structures Table

There is a Related Structures table for each feature in the Expected Features table and each compound in the Expected Compounds per File table. The Related Structures table shows the structure of the product compound generated by the dealkylation reaction.

* To open the Related Structures table

- 1. Open a result file from a targeted analysis (Find Expected Compounds node).
- 2. Select a row in one of these tables-Expected Features or Expected Compounds per File.

- 3. Click Show Related Tables.
- 4. Click the **Related Structures** tab.

Table 113 describes the columns in the Related Structures table.

Table 113. Related Structures table

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Parent Compound	Displays the user-specified parent compound or compounds.
Formula	Displays the chemical formula of the parent compound or the theoretical transformed compounds.
Molecular Weight	Displays the molecular weight (MW) of the parent compound or theoretical reaction product.
Dealkylated	Displays an X if the parent compound has undergone a dealkylation reaction.
Composition Change	Displays the composition change caused by any dealkylation reaction.
Structure	If a compound is the result of transformations that include a dealkylation step, this column displays the product of the dealkylation step. Otherwise, this column displays the structure of the parent compound.

Transformations Table

Use the Transformations table to review the transformations for each formula in the Expected Formulas table. Figure 7 on page 23 shows the relationship between the Expected Formulas table and the Transformations table.

Table 114 describes the columns in the Transformations table.

Table 114. Transformations table (Sheet 1 of 2)

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Name	Displays the name of the transformation.
Phase	Displays the Phase assignment for the transformation.
Leaving Group	Displays the leaving group for the transformation.
Arriving Group	Displays the arriving group for the transformation.

Column	Description
ΔMass [Da]	Displays the neutral mass shift for the transformation in daltons.
Order	Displays when the transformation was applied in the reaction pathway.
	Range: An integer from 1 to the user-specified maximum number of steps.

Table 114. Transformations table (Sheet 2 of 2)

Compound Detection Result Tables

For information about the result tables for the Compound Detection nodes, see these topics, listed in alphabetical order:

- Compounds Table
- Compounds per File Table
- Features Table
- Filled Gaps Table
- Labeled Features Table
- Labeled Compounds per File Table
- Similar Compounds (Beta) Related Table

Compounds Table

Use the Compounds table to review the unknown compounds found across the input file set.

The Group Compounds Node creates the Compounds table.

Clicking a row in the Compounds table displays XIC traces for the selected unknown compound, with one XIC trace for each input file. Each XIC trace is the summation of all the related ion traces. The summed XIC trace is made up of the data points with the highest intensity at each time point.

Tip To view the related traces for each XIC trace, do the following:

- 1. Select a row in the Compounds table.
- 2. In the first set of related tables, select a row in the Compounds per File table.
- 3. In the second set of related tables, open the Features table.

For information about exporting the compounds as a TXT or CSV file, see "Exporting the Tabular Results to an External File" on page 277. To edit the annotations, see "Editing Compound Annotations" on page 259.

Table 115 describes the columns in the Compounds table, with the additional columns in the related Compounds tables described at the end.

Column	Description
Checked	Specifies the rows that you want to display in the result table and in reports after you apply result filters.
Structure (hidden)	Displays the structure of the compound.
	The structure field is populated if any of the searches returns a structure or you edit the annotations.
Name	When the processing workflow includes a search node, this column displays the compound name from the best match in the searched databases.
Formula	Displays the predicted chemical formula for the neutral compound. For the analysis to predict and display a chemical formula for all compounds in the table, the processing workflow must include the Predict Compositions Node and the Assign Compound Annotations Node. When the processing workflow includes any of the search nodes, the Assign Compound Annotations node assigns the formula by using the specified priority for the data sources. The default priority for the data sources is (1) mzCloud Search, (2) Predicted Compositions, (3) Mass List Match, (4) ChemSpider Search.

Table 115. Compounds table (Sheet 1 of 8)

Column	Description
Annotation Source	Displays colorized rectangles that represent the match status for the selected compound from the search nodes in the processing workflow. The expanded column heading displays the annotation sources.
	The Assign Compound Annotations node determines the validity of the annotations from the Predict Compositions node and the search nodes in the processing workflow.
	Possible states for each annotation source:
	(■) Green—Full Match—The current formula and structure annotations match the best available item from the particular source (online database or local mass list).
	(□)Gray—No Results—Retrieved no data from the particular source.
	(D) Orange—Not the Top Hit—Current compound annotation matches one of the hits, but not the top one.
	() Orange—Partial Match—Only the formula for the current compound annotation matches the items retrieved from the particular source.
	(D) Orange—Unused—Retrieved items from the particular source, but did not assign any annotations.
	() Red—Invalid mass—The best available item from the particular source has a molecular weight that does not match the molecular weight of the compound within the specified mass tolerance.
	() Red—No match—The particular source does not have an item that matches the current annotations for the compound.
FISh Coverage	Displays the FISh Coverage score for a custom annotation (see "Editing Compound Annotations" on page 259 and "FISh Scoring for Targeted Compounds and Proposed Structures" on page 36).
Molecular Weight	Displays the molecular weight of the neutral compound.
RT [min]	Displays the retention time of the chromatographic peak for the compound.
RT Tolerance [min] (hidden)	Displays the retention time tolerance specified in the Group Compounds node.

 Table 115. Compounds table (Sheet 2 of 8)

Table 115. Compounds table (Sheet 3 of 8)

Column	Description
Area (Max.)	Displays the maximum chromatographic peak area from all of the input files for compounds with the same retention time and molecular weight (within the user-specified RT and mass tolerances).

The Normalize Areas Node and an input file set with QC samples add these columns to the result table: #Usable QC, RSD QC Areas [%], RSD QC Areas [%], QC Fill Status, and Norm. Areas.

For information about the batch normalization process, see "Quality Control Samples for Batch Normalization" on page 34.

To view the area correction for a compound, select the compound and choose **View** > **Compound Area Correction** (see "Viewing the QC-Based Compound Area Correction Curves" on page 367).

# Usable QC	Displays the number of usable QC samples (see Quality Control Samples for Batch Normalization).
RSD QC Areas [%]	Displays the relative standard deviation of the peak areas for the compound across the QC samples before area correction.
RSD Corr. QC Areas [%]	Displays the relative standard deviation of the peak areas for the compound across the QC samples after area correction.
QC Fill Status (hidden)	Displays a status rectangle for each QC sample. Possible states:
	Green—Filled by re-detected peak
	Gray—N/A
	Orange—Filled by matching ion
	Blue—Filled by simulated peak
Norm. Areas	(Normalize Areas node) Displays the normalized peak areas for the compound in each input file.
#ChemSpider Results	(Search ChemSpider Node) Displays the number of matching compounds found by the ChemSpider search for the current composition or molecular weight.
	Use the related ChemSpider Results Table to investigate the matching compounds.
Mass Defect	(Calculate Mass Defect node) Displays the compound's calculated mass defect.

Column	Description
#mzCloud Results	(Search mzCloud Node) Displays the number of matching compounds found by the mzCloud search for this molecular weight.
	Use the related mzCloud Results Table to investigate the matching compounds.
#Similarity Results (hidden)	Displays the number of similarity results from the Apply mzLogic node.
#mzVault Results	(Search mzVault Node) Displays the number of matching compounds found by the mzVault search for this molecular weight.
	Use the related mzCloud Results Table to investigate the matching compounds.
mzCloud Best Match	(Search mzCloud Node) Displays the best match score from the mzCloud identity search for the compound.
	Use the related mzCloud Results Table to investigate the matching compounds.
mzCloud Best Sim Match	(Search mzCloud Node) Displays the best similarity score from the mzCloud similarity search for the compound.
	Available when the Similarity Search parameter for the Search mzCloud node is set to Similarity Forward or Similarity Reverse.
	With the factory default setting, this column is the rightmost column in the table. Applying the Identification layout moves thi column to the right of the mzCloud Best Match column.
	Use the related mzCloud Results Table to investigate the matching compounds.
mzCloud Library Matches	Displays colorized rectangles that represent the match status for the compound in the selected mzCloud databases.
	Possible states for each database:
	() Green—Single match found
	(□)Gray—No matches found
	() Red—Multiple matches found

 Table 115. Compounds table (Sheet 4 of 8)

Column	Description
#Pathways (for any of the Map to Pathway nodes)	Displays the number of pathways that include the current compound.
	See Map to Metabolika Pathways Node, Map to BioCyc Pathways Node, or Map to KEGG Pathways Node.
Pathways	Displays whether the current compound is present in the named pathway.
	To display the pathway names, click the expand icon to the right of the column name.
	See Map to Metabolika Pathways Node, Map to BioCyc Pathways Node, or Map to KEGG Pathways Node.
Mass List Matches	(Search Mass Lists Node) Indicates the match status for each mass list.
	() Green—Single match found
	() Red—Multiple matches found
	() Gray—No matches found
	Use the related Mass List Search Results Table to investigate the matching compounds.
mzVault Library Matches	(Search Mass Lists Node) Indicates the match status for each library.
	() Green—Single match found
	() Red—Multiple matches found
	(□) Gray—No matches found
	Use the related mzVault Results Table to investigate the matching compounds.
Class Coverage	Displays the class coverage score for the compound. Expand the header to display the names of the compound class libraries.
#Adducts (hidden)	Displays the number of adduct ions that the analysis found for the compound.
Pattern Matches	(Pattern Scoring Node) Displays an orange rectangle when the compound matches the specified isotope pattern.
	Use the related Matched Patterns Table to investigate the matched isotope pattern.

 Table 115. Compounds table (Sheet 5 of 8)

Column	Description	
Labeling Status (per file)	The Analyze Labeled Compounds Node evaluates the measured isotope pattern versus the fitted isotope pattern (for the expected isotopologues) to determine the presence of contaminating masses. It also evaluates the distribution of the measured exchang rates for the expected isotopologues. If the distribution is not continuous, for example, if the compound has three exchangeab atoms and chromatographic peak area for the M+2 isotopologue significantly less than the chromatographic peak area for the M+ and M+3 isotopologues, the node assigns an Irregular Exchange status to the input file.	
	These flags indicate the following states:	
	() Red—Contaminating Mass—The average exchange for the unlabeled sample is above the 0.1 threshold.	
	() Orange—Low Pattern Fit—The measured pattern significantly differs from the fitted pattern. The SFit value is below the threshold of 20%, the Fitted Coverage value is below the threshold of 60%, or the Measured Coverage value is below threshold of 60%. To review these values, see the Labeled Feature table.	
	() Blue—Irregular Exchange—The isotopologue exchange rate are discontinuous; for example, there is a significant valley in the exchange rates profile. This might indicate an incorrect analysis of a special type of kinetics. However, if this is the typical behavior expected for your experiments, consider changing the setting for Mark Irregular Exchanges in the Analyze Labeled Compounds node to False.	
	Irregular Distribution 100 100 100 100 100 100 100 10	
	(■) Green—No Warnings—The measured isotope patterns and the exchange rates are within acceptable limits.	
	(\square) Gray—Compound was not detected in this sample.	
Avg. Exchange	Average number of atoms exchanged for compound detected in input file.	

 Table 115. Compounds table (Sheet 6 of 8)

Column	Description
Rel. Exchange [%]	Average exchange relative to the maximum exchange rate.
	100 × Average Exchange/Max. Exchange
MS2	Displays whether the analysis found data-dependent fragmentation scans for the compound.
	(■) Red—No MSn—There are no available MSn scans.
	(■) Green—ddMS2 for preferred ion—There is at least one data-dependent MS2 scan for the preferred adduct ion.
	(■) Blue—ddMS2 for other ion—There is at least one data-dependent MS2 scan, but the scans are not for the preferred adduct ion.
	(D)Orange—DIA only—Only data independent [all ions fragmentation (AIF)] scans are available.
Background	(Mark Background Compounds Node–Unknown Compounds) Displays a selected or clear check box that indicates whether the compound was also found in the Blank sample above the user-specified Sample/Blank or Blank/Sample level.
	 Selected—Indicates that the compound is a background compound.
	• Clear—Indicates that the compound is not a background compound.
Area (one column for	Displays an area column for each input file.
each input file) (hidden)	To display the areas for each input file, click the expand icon to the right of the column name.
Gap Status (hidden)	(Fill Gaps node)
	(■) Purple—Indicates some missing ions (but not a full gap).
	(□) Gray—Indicates a full gap.
	(■) Green—Indicates a compound without any gap.

Table 115. Compounds table (Sheet 7 of 8)

Table 115. Compounds table (Sheet 8 of 8)

Column Description	
	e following topics for information about the differential analysis columns and escriptive statistics columns.

- See "Descriptive Statistics Columns" on page 435 for information about the following columns: Min. Area, Median Area, Mean Area, Q1 Area, Q3 Area, Area CV%, and Area SD.
- See "Differential Analysis Columns" on page 432 for information about the following columns: Group Areas, Ratio, and Log2 Fold Change.

The following three columns appear in the related Compounds table for the main KEGG Pathways table. Each column contains an ordered list for the same KEGG compounds— that is, the first item in the KEGG Compound IDs column corresponds to the first item in the KEGG Compound Names column and the first item in the KEGG Compound Formula column, and so on.

KEGG Compound IDs	Displays a list of the KEGG compound IDs in ascending order from left to right.
KEGG Compound Names	Displays a list of the KEGG compound names.
KEGG Compound Formulas	Displays a list of the KEGG compound formulas.
The following two colur Pathways table and the s	nns appear in the related Compounds table for the main KEGG search results tables.
Max. ∆Mass [Da]	Displays the maximum mass difference between the measured mass for the compound in the Compounds table and the theoretical mass of the compound, in daltons.
Max. ΔMass [ppm]	Displays the maximum mass difference between the measured mass for the compound in the Compounds table and the theoretical mass of the compound, in parts per million.
This column appears in table.	the Compounds table that is related to the mzCloud Search Result
Scan Number	Displays the scan number of the scan that matches the reference

scan in the mzCloud database.

Compounds per File Table

Use the Compounds per File table to review the compounds detected in each input file.

Table 116 describes the columns in the Compounds per File table.

Table 116.Compounds per File table

Column	Description	
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.	
	Clear: False Selected: True	
Molecular Weight	Displays the molecular weight of the neutral compound.	
RT [min]	Displays the retention time at the chromatographic peak apex for the adduct ion that contributes the most area to the chromatographic peak.	
FWHM [min]	Displays the width of the chromatographic peak at its half-height. Use this value to determine the best RT tolerance for peak grouping.	
Max #MI	Displays the number of matching isotope peaks.	
#Adducts	Displays the number of adduct ions.	
Area	Displays the chromatographic peak area in counts * minutes.	

Features Table

Use the Features table to review the features detected across the input file set.

The Detect Compounds Node adds the Features table to the result file.

Table 117 describes the columns in the Features table.

Table 117. Features table (Sheet 1 of 2)

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Ion	Displays the ion definition of the molecular ion adduct.
Charge	Displays the charge on the ion.
Molecular Weight	Displays the molecular weight of the unknown compound.
m/z	Displays the mass-to-charge ratio of the ion.
RT [min]	Displays the retention time in minutes of the chromatographic peak that contains the unknown compound ion.

Column	Description
#MI	Displays the number of matching isotopes for the unknown compound ion.
Intensity (hidden)	Displays the intensity of the ion.
Area	Displays the area of the chromatographic peak that contains the unknown compound ion.
Parent Area [%]	Displays the chromatographic peak area of the current peak as a percentage of the total chromatographic peak area for the parent compound (which is the compound selected in the Compounds per File table) per input file.
File ID	Displays the integer that the application assigned to the input file.
Study File ID	Displays the study file ID of the input file.

Table 117	. Features table	(Sheet 2 of 2)
-----------	------------------	----------------

Filled Gaps Table

Use the Filled Gaps table to review the chromatographic peaks (Full Gap or Missing Ion) that the Fill Gaps node finds. The Filled Gaps table is related to the Compounds table.

By default, the Filled Gaps table is hidden. To display the table, click the **Select Table Visibility** icon, ^(III), select the **Filled Gaps** check box, and click **OK**.

The Gap Status column, hidden by default in the Compounds table, identifies the chromatographic peaks that the Fill Gaps node finds as Full Gap or Missing Ion peaks and those that the standard peak detection process finds as No Gap peaks.

Table 118 describes the columns in the Filled Gaps table.

Table 118. Filled Gaps table (Sheet 1 of 2)

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Ion Description	Displays a description of the adduct where M is the neutral compound.
Charge	Displays the ion's charge.
Exp. m/z	Displays the <i>m/z</i> value for the adduct ion found in the other input files. This value is based on the adduct type and the compound's molecular weight.
Exp. RT [min]	Displays the average retention time of the adduct ion found in the other input files.

Column	Description
Exp. Left RT [min]	Displays the average peak start time of the detected chromatographic peaks for the adduct ion in all the input files included in the analysis.
Exp. Right RT [min]	Displays the average peak end time of the detected chromatographic peaks for the adduct ion in all the input files included in the analysis.
Exp. FWHM [min]	Displays the average peak width at the peak's half height (full width at half maximum) for the adduct ion in all the input files included in the analysis.
Area	Displays the area of the chromatographic peak found by the Fill Gaps node.
Fill Status	The possible states for the chromatographic peak detected or redrawn by the Fill Gaps node are as follows:
	• () Green (Filled by Simulated Peak or Filled by Re-detected Peak)—The chromatographic peak was detected with the PPD algorithm (set to a lower threshold than in the Detect Compounds node) or the peak was simulated with a Gaussian fit algorithm.
	• () Blue (Filled by Spectrum Noise)—The gap was replaced with a chromatographic peak based on the spectrum noise level.
	• () Orange (Filled by Matching Ion)—The gap was replaced with a chromatographic peak for a matching ion.
File ID	Displays the integer that the application assigned to the input file.
Study File ID	Displays the study file ID of the input file

Table 118. Filled Gaps table (Sheet 2 of 2)

Labeled Features Table

Table 119 describes the columns in the Labeled Features table. The labeled features are the labeled adduct ions.

Table 119. Labeled Features table (Sheet 1 of 2)

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Ion	Displays the ion definition of the molecular ion adduct.
Charge	Displays the charge on the ion.

Description	
Displays the molecular weight of the unknown compound.	
Displays the mass-to-charge ratio of the ion.	
Displays the retention time in minutes of the chromatographic peak that contains the unknown compound ion.	
Displays the width of the chromatographic peak for the adduct ion at its half-height in minutes.	
Displays the intensity of the ion.	
Displays the area of the chromatographic peak that contains the unknown compound ion.	
Displays the chromatographic peak area of the current peak as a percentage of the total chromatographic peak area for the parent compound (compound selected in the Compounds per File table) per input file.	
Displays the maximum number of atoms considered for isotopologue evaluation.	
Displays the average number of atoms exchanged for the ion.	
Displays the relative number of atoms exchanged versus the maximum number of exchangeable atoms for the feature's elemental composition.	
Displays the exchange rate for individual isotopologues.	
Displays the number of matching isotopes for the unknown compound ion.	
Displays the spectral similarity score between the measured and theoretical isotope pattern.	
Displays the how well the intensities of the fitted isotope pattern match those of the theoretical isotope pattern.	
Displays the how well the intensities of the measured isotope pattern match those of the theoretical isotope pattern.	
Displays the integer that the application assigned to the input file.	
Displays the study file ID of the input file (see "Removing Input Files or Updating Their Location" on page 105).	

 Table 119.
 Labeled Features table (Sheet 2 of 2)

Labeled Compounds per File Table

Table 120 describes the columns in the Labeled Compounds per File table. The related table displays details about the selected compound in the higher level table. To view the compound name for a component in the main Labeled Compound per File table, open its related Compounds table.

Column	Description	
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.	
Molecular Weight	Displays the molecular weight of the unknown compound.	
RT [min]	Displays the retention time in minutes of the chromatographic peak that contains the unknown compound ion.	
FWHM	Displays the width of the chromatographic peak for the adduct ion at its half-height in minutes.	
Max. #MI	Displays the maximum number of matching isotope peaks.	
#Adducts	Displays the number of adducts (features).	
Area	Displays the area of the chromatographic peak that contains the unknown compound ion.	
Max. Exchange	Shows the maximum number of exchangeable atoms that the analysis considered.	
Avg. Exchange	Displays the average number of atoms exchanged for the ion.	
Rel. Exchange [%]	Displays the relative number of atoms exchanged versus the maximum number of exchangeable atoms.	
Status	The Analyze Labeled Compounds Node evaluates the measured isotope pattern versus the fitted isotope pattern (for the expected isotopologues) to determine the presence of contaminating masses. It also evaluates the distribution of the measured exchange rates for the expected isotopologues.	
	For information about the status flags, see "Labeling Status" on page 407.	
Exchange Rate [%]	Displays the contribution of individual isotopologues to the final measured pattern.	
File ID	Displays the integer that the application assigned to the input file.	
Study File ID	Displays the study file ID of the input file.	

Similar Compounds (Beta) Related Table

Use the Similar Compounds (Beta) table to review the compounds that the application connected to the detected compound in the Compounds table.

The Similar Compounds (Beta) table is generated by the Generate Molecular Networks node.

Table 121 describes the columns in the Similar Compounds table.

Table 121. Similar Compounds table (Sheet 1 of 2)

Column	Description		
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.		
Direction	Displays the direction in which the transformation was applied, and defines which compound is a substrate and which is a product.		
	Forward—The selected compound acts as a substrate to which the transformation was applied. The connected similar compound is a product of the transformation.		
	Reverse—The selected compound is a product of the transformation. The connected similar compound acts as a substrate to which the transformation was applied.		
Mass Shift [Da]	Displays the difference between the measured molecular weights of the two compounds.		
Composition Change	Displays the difference between the elemental compositions of the two compounds.		
Transformations	Displays the name of assigned transformation pathway—that is, this column lists the names of the individual steps. Displays "Isomer" when the elemental composition of the substrate and product compounds are identical.		
Transformation Mass [Da]	Displays the total theoretical mass of the assigned transformation steps in Daltons.		
# Fragments	Displays the number of fragments available for the selected compound.		
MSn Score	Displays the final spectral similarity score between the two compounds as the average of the forward and reverse coverages.		
Forward Cov. [%]	Displays the relative number of matched centroids in the product compound's MSn spectra.		
Forward Matches	Displays the number of matched centroids in the product compound's MSn spectra.		

Column	Description
Reverse Cov. [%]	Displays the relative number of matched centroids in the substrate compound's MSn spectra.
Reverse Matches	Number of matched centroids in the substrate compound's MSn spectra.

Table 121. Similar Compounds table (Sheet 2 of 2)

Compound Identification Result Tables

For information about the result tables for the Compound Identification workflow nodes, see these topics:

- ChemSpider Results Table
- Mass List Search Results Table
- mzCloud Results Table
- mzVault Results Table
- Predicted Compositions Table

ChemSpider Results Table

Use the ChemSpider Results table to review the compounds found in the ChemSpider databases. The Search ChemSpider Node creates the ChemSpider Results table.

* To open the ChemSpider information for a specific ChemSpider hit

Click the link in the CSID column.

The ChemSpider web page for the selected hit opens in your default web browser.

Table 122 describes the columns in the ChemSpider Results table.

Table 122.	ChemSpider	Results table	(Sheet 1 of 2)
------------	------------	---------------	----------------

Column	Description	
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.	
Name	Displays the name of the compound hit in the ChemSpider database.	
Structure	Displays the molecular structure of the compound.	
Formula	Displays the chemical formula of the compound.	

Column	Description	
Molecular Weight	Displays the molecular weight of the compound to five decimal places.	
CSID	Displays the ChemSpider identification number.	
#References	Displays the number of references for the compound in the ChemSpider database.	
Additional hidden colum	ins	
SMILES (hidden)	Displays the compound's molecular structure by using short ASCII strings. SMILES stands for simplified molecular input line entry System.	
InChi (hidden)	Displays the international chemical identifier for the compound.	
#Data Sources (hidden)	Displays the number of ChemSpider data sources that include the compound.	
#PubMed References (hidden)	Displays the number of PubMed references for the compound. You can use a PubMed reference to access the scientific literature	
AlogP (hidden)	Displays the AlogP value for the compound. The AlogP value is a measure of the compound's hydrophobicity, where P is the partition coefficient.	
#RSC (hidden)	Displays the number of Royal Society of Chemistry references fo the compound.	
XlogP (hidden)	Displays the XlogP value for the compound. The XlogP value is a measure of the compounds hydrophobicity, where P is the partition coefficient.	
Additional columns in th	e related ChemSpider Results table	
Compound Match	Indicates the match status between the current item and the assigned compound annotation. For color-coding information, see "Annotation Source" on page 403.	
ΔMass [Da]	Displays the mass difference in daltons between the search mass and the mass of the matching compound in the ChemSpider database.	
ΔMass [ppm]	Displays the mass difference in ppm between the search mass and the mass of the matching compound in the ChemSpider database	
SFit[%]	Displays the spectral similarity score from the Apply Spectral Distance node.	

 Table 122.
 ChemSpider Results table (Sheet 2 of 2)

Mass List Search Results Table

Use the Mass List Search Results table to review the compounds in the selected mass lists that match the compounds detected by the Detect Compounds node.

The Search Mass Lists Node creates the Mass List Search Results table.

The main Mass List Search Results table displays all of the compounds in the selected mass lists that match the compounds detected by the untargeted analysis. The related Mass List Search Results table lists information about the compound selected in the main Compounds Table.

Table 123 describes the columns in the Mass List Search Results table.

Column	Description	
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.	
Formula	Displays the elemental composition of the compound in the mass list.	
Molecular Weight	Displays the molecular weight of the compound in the mass list.	
RT [min]	Displays the chromatographic retention time (if available) of the compound in the mass list.	
Structure	Displays the structure (if available) of the compound in the mass list.	
Name	Displays the name of the compound in the mass list.	
Annotation	Displays additional information about the compound in the mass list.	
Reference List Name	Displays the name of the mass list that contains the matching compound.	
Additional columns in the related Mass List Search Results table		
Compound Match	Indicates the match status between the current item and the assigned compound annotation. For color-coding information, see "Annotation Source" on page 403.	
Max. ΔMass [Da]	Displays the maximum mass difference between the measured mass for the compound in the Compounds table and the theoretical mass of the compound, in daltons.	

 Table 123. Mass List Search Results table (Sheet 1 of 2)

Column	Description
Max. ΔMass [ppm]	Displays the maximum mass difference between the measured mass for the compound in the Compounds table and the theoretical mass of the compound, in parts per million.
SFit[%]	Displays the spectral similarity score from the Apply Spectral Distance node.

Table 123. Mass List Search Results table (Sheet 2 of 2)

mzCloud Results Table

Use the mzCloud Results table to view the results of an mzCloud search.

The Search mzCloud Node creates the mzCloud Results table.

Selecting a row in the mzCloud Results table displays a mirror plot with the selected fragmentation scan on the top and the matched reference scan from the mzCloud database on the bottom. Figure 121 shows a mirror plot with structure annotations. The centroids for the matching fragments are displayed as green sticks with a green circle at the end. Red circles on the *x*-axis indicate the m/z values of the missing fragments.

Note By default, the Search mzCloud node does not annotate the matching fragments. To annotate the matching fragments, you must enable this feature—that is, you must set the node's Annotate Matching Fragments parameter to True.

Figure 121. Mass Spectrum view with an annotated mirror plot

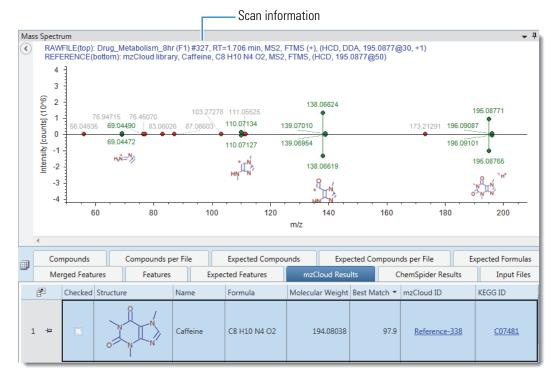


Table 124 describes the columns in the mzCloud Results table.

Table 124	. mzCloud	Results	table	(Sheet 1	of 2)
-----------	-----------	---------	-------	----------	-------

Column	Description	
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.	
Structure	Displays the structure of the matching compound.	
Name	Displays the compound name in the mzCloud database.	
Formula	Displays the elemental composition formula of the matching compound.	
Molecular Weight	Displays the molecular weight of the matching compound.	
Best Match	Displays the match value (0–100%) between the best fragmentation scan for a compound (across the input files) and the matching mzCloud spectrum (mzCloud ID). The related Compounds tables lists the scan number of the best fragmentation scan.	
Best Sim. Match	Displays the best match score for a library spectrum to a fragmentation scan across the input files. The related Compounds tables lists the scan number of the similar fragmentation scan.	
mzCloud ID	Displays the mzCloud ID (<i>Database-Number</i>) for the matching compound. Clicking this link opens the mzCloud database to the matching compound's reference spectrum.	
KEGG ID	Displays the KEGG ID for the compound in the KEGG database.	
Compound Class	Displays the mzCloud compound classes that include the compound.	
mzCloud Library	Displays the name of the database where the search found the matching spectrum.	
Additional columns in th	e related mzCloud Results table	
Compound Match	Indicates the match status between the current item and the assigned compound annotation. For color-coding information, see "Annotation Source" on page 403.	
Searched with ∆Mass [Da]	Displays the maximal difference between the theoretical and measured mass in Da.	
mzCloud Library	Displays the name of the mzCloud database where the search found the matching scan.	
Intensity Threshold	Displays the relative intensity threshold for the search spectrum.	
ΔMass [Da]	Displays the difference between the theoretical mass of the matching compound and the observed mass of the unknown compound in daltons.	

Column	Description	
ΔMass [ppm]	Displays the difference between the theoretical mass of the matching compound and the observed mass of the unknown compound in parts per million.	
Туре	Displays the search type: Identify or Similarity.	
Scan Number	Displays the scan number from the input files that contains the best matching fragmentation spectrum.	
Match	Displays the match value (0–100%) between the specified scan number from the input files and the matching spectrum in the mzCloud database.	

Table 124. mzCloud Results table (Sheet 2 of 2)

Related Topics

• Search mzCloud Node

mzVault Results Table

Use the mzVault Results table to review the results of an mzVault search.

The Search mzVault Node creates the mzVault table.

Table 125 describes the columns in the mzVault Results table.

Table 125. mzVault Results table (Sheet 1 of 2)

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Structure	Displays the structure of the matching compound.
# Related Compounds	Displays the number of unknown compounds with the same putative elemental composition but different retention times.
mzVault ID	Displays the ID number of the matching compound in the mzVault library.
Name	Displays the name of the matching compound.
Formula	Displays the elemental composition formula of the matching compound.
Molecular Weight	Displays the molecular weight of the matching compound.
Best Match	Displays the match value (0–100%) between the best fragmentation scan for a compound (across the input files) and the matching mzVault spectrum. The related Compounds table lists the scan number of the best fragmentation scan.

Column	Description
mzVault Library	Displays the name of the mzVault library where the analysis found a matching compound.
ChemSpider ID	Displays the ChemSpider ID number for the matching compound. Clicking this link opens the ChemSpider database to the compound's record.
Compound Class	Displays the mzVault compound class that includes the compound.
mzCloud ID	Displays the mzCloud ID for the matching compound in the mzCloud database. Clicking this link opens the mzCloud database to the compound's record.
KEGG ID	Displays the KEGG ID for the matching compound in the KEGG database. Clicking this link opens the KEGG database to the compound's record.
Additional columns in t	he related mzVault Results table
Compound Match	Indicates the match status between the current item and the assigned compound annotation. For color-coding information, see "Annotation Source" on page 403.
ΔMass [Da]	Displays the difference between the theoretical mass of the matching compound and the observed mass of the unknown compound in daltons.
ΔMass [ppm]	Displays the difference between the theoretical mass of the matching compound and the observed mass of the unknown compound in parts per million.
Scan Number	Displays the scan number from the input files that contains the best matching spectrum.
Match	Displays the match value (0–100%) between the specified scan number from the input files and the matching spectrum in the mzVault database.

Table 125. mzVault Results table (Sheet 2 of 2)

Predicted Compositions Table

Each compound in the Compounds table has a related Predicted Compositions table.

Use the Predicted Compositions table to review the possible chemical formulas for the selected compound in the Compounds table. The Predicted Compositions tables lists the possible chemical formulas based on the compound's molecular weight.

The Predict Compositions Node creates the Predicted Compositions table.

Table 126 describes the columns in the Predicted Compositions result table.

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Compound Match	Displays whether the predicted composition matches the currently assigned compound annotation.
	(Red rectangle—No Match
	() Green rectangle—Full Match
Formula	Displays the predicted elemental composition.
Molecular Weight	Displays the molecular weight.
ΔMass [Da]	Displays the difference between the theoretical mass and the measured mass in daltons.
ΔMass [ppm]	Displays the difference between the theoretical mass and the measured mass in ppm.
RDBE	Displays the rings and double bonds equivalent value for the predicted composition.
H/C	Displays the ratio of hydrogen to carbon atoms in the predicted composition.
Rank	Displays the rank order of each composition.
#Matched Iso.	Displays the number of matching isotopes.
#Missed Iso.	Displays the number of isotopes that were missing in the measured isotope pattern as compared to the theoretical pattern for the predicted composition.
#Matched Fragments	When the Use Fragments Matching algorithm is turned on in the Predict Compositions node, this column displays the number of centroids (m/z values) in the best MS2 scan that match possible fragments (mass values from a subset of the elemental compositions in the predicted composition).
SFit [%]	Displays the spectral similarity score between the theoretical and the measured isotope pattern as a percentage.
	The SFit [%] score = $(1-SD) \times 100$
SD (hidden)	Displays the spectral distance score. A lower SD score corresponds to a higher SFit [%] score.
	Range: 0 to 1

Table 126. Predicted Compositions table (Sheet 1 of 2)

Column	Description
Pattern Cov. (%)	Displays the summed intensity of the matching isotope peaks in the measured MS1 spectrum relative to the summed intensity of the theoretical isotope pattern.
	Summed intensity of the matching isotope peaks \times 100 Summed intensity of the theoretical isotope pattern
	Provides a quantitative measure of how well the measured isotope pattern matches the theoretical isotope pattern.
	Note Because the base peak (leftmost peak) is typically responsible for most of the pattern intensity, even a small decrease in the percent coverage might be important. For example, a missing peak for an isotope with two ¹³ C atoms might cause only a small decrease in the summed intensity of the measured isotope pattern.
MS Cov. (%)	Displays the summed intensity of matching isotope peaks in the measured pattern relative to the summed intensity of all the peak in the measured pattern.
	$\frac{\text{Summed intensity of the matching isotope peaks} \times 100}{\text{Summed intensity of all the peaks in the measured pattern}}$
	IMPORTANT Low values for all of the candidates might indicate an overlapping pattern rather than a lack of good matches.
MSMS Cov. (%)	Displays the summed intensity of the matched fragment peaks relative to the summed intensity of all the fragment peaks in the selected fragmentation scan.
	Note Low values for all of the candidates might indicate a contaminating compound within the isolation window for the fragmentation scan.

Table 126. Predicted Compositions table (Sheet 2 of 2)

Pathway Mapping Result Tables

For information about the result tables for the Pathway Mapping nodes, see these topics:

- BioCyc Pathways Table
- BioCyc Results Table
- KEGG Pathways Table
- Metabolika Pathways Table
- Metabolika Results Table

BioCyc Pathways Table

Use the BioCyc Pathways table to review all the mapped pathways in the result file.

For more information, see "Viewing BioCyc Pathways" on page 362.

Table 129 describes the columns in the BioCyc Pathways main table. The BioCyc Pathways table lists all the BioCyc pathways that include at least one of the compounds detected by the untargeted search.

Table 127.BioCyc Pathways table

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Pathway ID	Displays the identification number of the mapped BioCyc pathway.
Pathway Name	Displays the name of the mapped BioCyc pathway that includes at least one compound in the Compounds table.
#Referenced Compounds	Displays the number of detected compounds that are referenced in the identified BioCyc pathway.
#Identified Compounds	Displays the number of compounds in the referenced pathway that the analysis detected in the input file set.
	To view the list of detected compounds, select a pathway and open the related BioCyc Results table.
Related BioCyc Pathways tabl	e only
BioCyc Compound IDs	Displays a list of the BioCyc compound IDs in ascending order from left to right.
BioCyc Compound Names (related table)	Displays a list of the BioCyc compound names.
BioCyc Compound Formulas (related table)	Displays a list of the BioCyc compound formulas.

BioCyc Results Table

Use the BioCyc Results table to review the compounds found in the mapped BioCyc database.

The Map to BioCyc Pathways Node creates the BioCyc Results table.

Table 128 describes the columns in the BioCyc Results table.

Table 128.	BioCyc Results table
------------	----------------------

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Structure	Displays the molecular structure of the compound.
Name	Displays the compound name in the BioCyc database.
Formula	Displays the chemical formula of the compound.
Molecular Weight	Displays the molecular weight of the compound to five decimal places.
Additional columns in t	he related BioCyc Results table for a compound
Compound Match	Indicates the match status between the current item and the assigned compound annotation. For color-coding information, see "Annotation Source" on page 403.
Max. ΔMass [Da]	Displays the maximum mass difference between the measured mass for the compound in the Compounds table and the theoretical mass of the compound, in daltons.
Max. ΔMass [ppm]	Displays the maximum mass difference between the measured mass for the compound in the Compounds table and the theoretical mass of the compound, in parts per million.
BioCyc ID	Displays the BioCyc identification number for the compound.
BioCyc DB	Displays the name of the BioCyc database that mapped the compound.
mzLogic Score	Displays the mzLogic score provided by the Apply mzLogic node.
SFit[%]	Displays the spectral similarity score from the Apply Spectral Distance node.

KEGG Pathways Table

Use the main KEGG Pathways table to review all of the mapped pathways that include compounds detected across the input file set.

The Map to KEGG Pathways Node creates the KEGG Pathways table.

Table 129 describes the columns in the KEGG Pathways table. The KEGG Pathways table lists all the KEGG pathways that include at least one of the compounds detected by the untargeted search.

Table 129.	KEGG	Pathways	table
------------	------	----------	-------

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Pathway ID	Displays the identification number of the mapped KEGG pathway.
Pathway Name	Displays the name of the mapped KEGG pathway that includes at least one compound in the Compounds table.
#Referenced Compounds	Displays the number of detected compounds that are referenced in the identified KEGG pathway.
#Identified Compounds	Displays the number of different KEGG compound IDs.
#mzCloud Results	Displays the number of compounds in the identified KEGG pathway that the mzCloud search identified.
KEGG Compound IDs (related table)	Displays a list of the KEGG compound IDs in ascending order from left to right.
KEGG Compound Names (related table)	Displays a list of the KEGG compound names.
KEGG Compound Formulas (related table)	Displays a list of the KEGG compound formulas.
Max. ΔMass [Da]	Displays the maximum mass difference between the measured mass for the compound in the Compounds table and the theoretical mass of the compound, in daltons.
Max. ΔMass [ppm]	Displays the maximum mass difference between the measured mass for the compound in the Compounds table and the theoretical mass of the compound, in parts per million.

Metabolika Pathways Table

Use the main Metabolika Pathways table to review all of the mapped pathways that include compounds detected across the input file set. Use the related Metabolika Pathways tables to review the results for specific compounds.

The Map to Metabolika Pathways Node creates the Metabolika Pathways table.

Table 130 describes the columns in the Metabolika Pathways table. The main Metabolika Pathways table lists all the Metabolika pathways that include at least one of the compounds detected by the untargeted search. The related table displays the search results for the selected compound in the Compounds table.

Table 130. Metabolika Pathways table

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Pathway Name	Displays the names of the mapped Metabolika pathways that include matching structures (by formula, mass, or both) for at least one compound in the Compounds table.
#Referenced Compounds	Displays the number of detected compounds that are referenced in the identified Metabolika pathway.
#Compounds in Pathway	Displays the total number of compounds in the pathway.
#Identified Compounds	Displays the number of different Metabolika compound IDs.
Metabolika Compound IDs (related table)	Displays a list of the Metabolika compound IDs in ascending order from left to right.
Metabolika Compound Names (related table)	Displays a list of the Metabolika compound names.
Metabolika Compound Formula (related table)	Displays the Metabolika compound formula.

Metabolika Results Table

Use the main Metabolika Results table to review the compounds found in the mapped Metabolika pathways. Use the related Metabolika results table to review the results for a specific compound.

The Map to Metabolika Pathways Node creates the Metabolika Results table.

Table 128 describes the columns in the Metabolika Results table.

 Table 131.
 Metabolika Results table (Sheet 1 of 2)

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Structure	Displays the molecular structure of the compound.
Name	Displays the compound name in the Metabolika database.
Formula	Displays the chemical formula of the compound.

Column	Description
Molecular Weight	Displays the molecular weight of the compound to five decimal places.
Additional columns in th	e Metabolika Results related table
Compound Match	Displays whether the Metabolika pathway structure is a full or partial match for the selected compound.
	()—Full Match
	(])—Partial Match
mzLogic Score	Displays the normalized score from the Apply mzLogic node for the structure based on the mzCloud similarity matches (Apply mzLogic node).
Original mzLogic Score	Displays the score from the Apply mzLogic node for the structure based on the mzCloud similarity matches
Max. ΔMass [Da]	Displays the maximum mass difference between the measured mass for the compound in the Compounds table and the theoretical mass of the compound, in daltons.
Max. ΔMass [ppm]	Displays the maximum mass difference between the measured mass for the compound in the Compounds table and the theoretical mass of the compound, in parts per million.
SFit[%]	Displays the spectral similarity score from the Apply Spectral Distance node.

Table 131. Metabolika Results table (Sheet 2 of 2)

Compound Scoring Tables

For information about the result tables for the unknown compound scoring nodes, see these topics:

- Compound Class Matches Table
- Matched Patterns Table

Compound Class Matches Table

The Compound Class Matches table is related to the Compounds table. Use the Compound Class Matches table to review the compounds classes that match the compound selected in the Compounds table.

The Compound Class Scoring Node adds the Class Coverage column to the Compounds table and creates the Compound Class Matches table, which is related to the Compounds table.

Table 132 describes the columns in the Compound Class Matches table.

Table 132.	Compound	Class	Matches table
------------	----------	-------	---------------

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Name	Displays the name of the library that contains the matching structure.
Description	Displays the user-specified description of the library.
FISh Coverage	Displays the FISh coverage score (see "FISh Scoring for Targeted Compounds and Proposed Structures" on page 36).
Class Coverage	Displays the number of matching centroids in the best fragmentation scan divided by the total number of fragments in the selected compound class libraries.
# Matched Fr.	Displays the number of library fragments that match the centroids in the best fragmentation scan for a compound.
# Missed Fr.	Displays the number of library fragments that do not match the centroids in the best fragmentation scan for a compound.

Figure 122 shows an annotated fragmentation scan of a detected compound. By comparing the m/z values of the centroids in the fragmentation scan against a compound class library with 11 structures, the application annotated 6 centroids in the fragmentation scan with matching structures from the library. The legend in the Mass Spectrum view lists the search library, the Class Coverage score, and the FISh Coverage score.

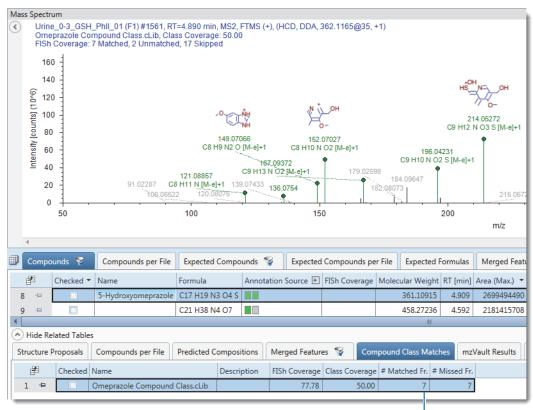


Figure 122. Fragmentation scan and compound class match result for a detected compound

#Matched fragments ----

Matched Patterns Table

The Matched Patterns table is related to the Compounds table. Use the Matched Patterns table to review how well the isotopic pattern matches the compound selected in the Compounds table.

The Pattern Scoring node adds the Matched Patterns table to the result file.

Table 133 describes the columns in the Matched Patterns table.

Table 133. Matched Patterns table (Sheet 1 of 2)

Column	Description
Checked	Use this column to select the rows that you want to display in the result table and in reports after you apply result filters.
Name	Displays the name or chemical formula of the compound.
SFit [%]	Displays the spectral fit for the isotope pattern to the chemical formula.

Description
Displays the spectral distance score.
Displays the summed intensity of the matching isotope peaks in the measured MS1 spectrum relative to the summed intensity of the theoretical isotope pattern.
$\frac{\text{Summed intensity of the matching isotope peaks} \times 100}{\text{Summed intensity of the theoretical isotope pattern}}$
Provides a quantitative measure of how well the measured isotope pattern matches the theoretical isotope pattern.
Note Because the base peak (leftmost peak) is typically responsible for most of the pattern intensity, even a small decrease in the percent coverage might be important. For example, a missing peak for an isotope with two ¹³ C atoms might cause only a small decrease in the summed intensity of the measured isotope pattern.
Displays the number of matching isotopes for the unknown compound.
Displays the integer that the application assigned to the input file.
Displays the study file ID of the input file (see "Removing Input Files or Updating Their Location" on page 105).

Table 133. Matched Patterns table (Sheet 2 of 2)

Differential Analysis Columns

When the analysis includes sample groups and group ratios and the processing workflow includes the Differential Analysis node, the following columns appear in the Compounds Table and the Expected Compounds Table: Group Areas, Group CV(%), Ratio, Log2 Fold Change, P-value, and Adj. P-value. In addition, the following columns appear in the Merged Features Table, Group Areas, Ratio, and Log2 Fold Change. The background colors of the table cells provide visual information about the numeric values in these columns.

Table 134 describes the columns from a differential analysis.

Table 134. Differential analysis columns (Sheet 1 of 3)

Column	Description
Group Areas	Displays the median chromatographic peak area for the compound in the sample group. To display the group names, click the expand icon to the right of the column heading.
	When the compound is not found, the cells have a gray background. The Differential Analysis node bins the group areas in the current result file on the Log10 scale, with one bin for each order of magnitude (1e4, 1e5, 1e6, and so on), and uses a different background color for each bin. The values in the lowest bin have a pale yellow background. The values in the highest bin have a green background.
	Color-coding:
	Lowest values:
	Highest values:
Group CV(%)	Displays the coefficient of variation for the groups. Groups with a hig degree of variation (20% or greater) have a red background.
	Color-coding:
	(20) Values equal to or greater than 20 have a red background.
	(5) Values from 0 to 19 have a white background.
Ratio	Displays the area ratio or ratios for the generated ratios.
	A ratio of 0 (0/X) has a dark purple background (0.000). An undefined ratio (X/0) (labeled as Infinity) has an orange background (Infinity).
	The cells for compounds with defined ratios greater than 0 have the following background colors:
	• Compounds with ratios between 0 and 0.5 are divided into five equal bins. The background color for the table cells is a progressively darker blue hue as the ratio approaches zero.
	• Compounds with ratios greater than 2 are divided into five equa bins. The background color for the tables cells is a progressively darker red hue as the ratio increases.
	0 . - 0
	0.5 2 0 < Ratio < 0.5 Ratio ≥ 2

Column	Description	
Log2 Fold Change	Displays the fold change (ratio) in the log base 2 scale.	
	A log2 fold change of Infinity has an orange background (<mark>Infinity</mark>). I the ratio is 0 (0/X) or –Infinity in the Log2 format, the background color is a dark purple (<mark>Infinity</mark>).	
	The cells for compounds with log2 fold change values between –infinity and infinity have the following background colors:	
	• Compounds with log2 fold change values more negative than – 1.00 are divided into five equal bins. The background color for the table cells is a progressively darker blue hue as the value becomes more negative.	
	• Compounds with log2 fold change values greater than 1.00 are divided into five equal bins. The background color for the table cells is a progressively darker red hue as the value increases.	
	Log2 Fold Change negative −1 1 positive ←	
P-value	Displays the p-value for the sample group calculated by running the Tukey HSD test (posthoc) after an analysis of variance (ANOVA) tes	
	The p-value is a number between 0 and 1.	
	Given the following hypotheses:	
	• Null hypothesis—There is no difference between the sample groups for the variable tested.	
	• Alternate hypothesis—There is a difference between the sample groups for the variable tested.	
	You can interpret the p-value as follows:	
	• A low p-value means that you can reject the null hypothesis with low probability of error that the alternate hypothesis is true.	
	• A high p-value means that you can accept the null hypothesis with a low probability of error that the alternate hypothesis is true.	

Table 134. Differential analysis columns (Sheet 2 of 3)

Column	Description
Adj. P-Value	(Differential Analysis node) Displays the adjusted p-value.
	The application adjusts p-values in cases of multiple testing. Multiple testing of a null hypothesis leads to higher probabilities of rejecting this null hypothesis by chance, and therefore the application corrects the whole set of hypotheses (for example, all detected compounds) as a function of the set size (for example, a set of 10 000 compounds has a stronger correction than one of only 1000). The application performs this correction by using the Benjamini-Hochberg algorithm for the false discovery rate.
P-value and Adj.	Color-binning for p-values and adjusted p-values:
P-value	• 1–0.05 [
	• 0.05–0.01 [] Orange
	• 0.01–0.005 [] Yellow
	• 0.005–0.001 [] Yellow green
	• <0.001 [] Green

Table 134. Differential analysis columns (Sheet 3 of 3)

Descriptive Statistics Columns

The Descriptive Statistics post-processing node adds the columns described in Table 135 to the Compounds Table and Expected Compounds Table. The descriptive statistics columns are hidden by default.

Column	Description
Min. Area	Displays the minimum peak area for the compound in the sample set.
Q1 Area	Displays the lower boundary of the first quartile (25%) area for the compound in the sample set.
Median Area	Displays the median area for the compound (MW×RT) in the sample set.
Q3 Area	Displays the upper boundary of the third quartile (75%) area for this compound in the sample set.
Mean Area	Displays the calculated average area for the compound (MW×RT) in the sample set.

Table 135. Descriptive statistics columns (Sheet 1 of 2)

Column	Description
Area SD	Displays the standard deviation of the peak areas for the compound (MW×RT) in the sample set.
Area CV [%]	Displays the coefficient of variation of the area for the compound (MW×RT) in the sample set.

Table 135. Descriptive statistics columns (Sheet 2 of 2)

11

Creating and Printing Reports

These topics describe how to create, preview, and print reports. You can use the standard report templates provided with the application or you can create your own custom report templates.

Contents

- Reporting Workflow
- Generating a Report with an Existing Report Template
- Creating a New Report Template
- Editing an Existing Report Template
- Editing Tools for the Report Template Page
- Section Report Items for a Report Template
- Property Settings for the Sections and Items in a Report Template
- Selecting the Paper Type, Print Width, Page Orientation, and Watermark for a Report Template
- Previewing and Printing a Report

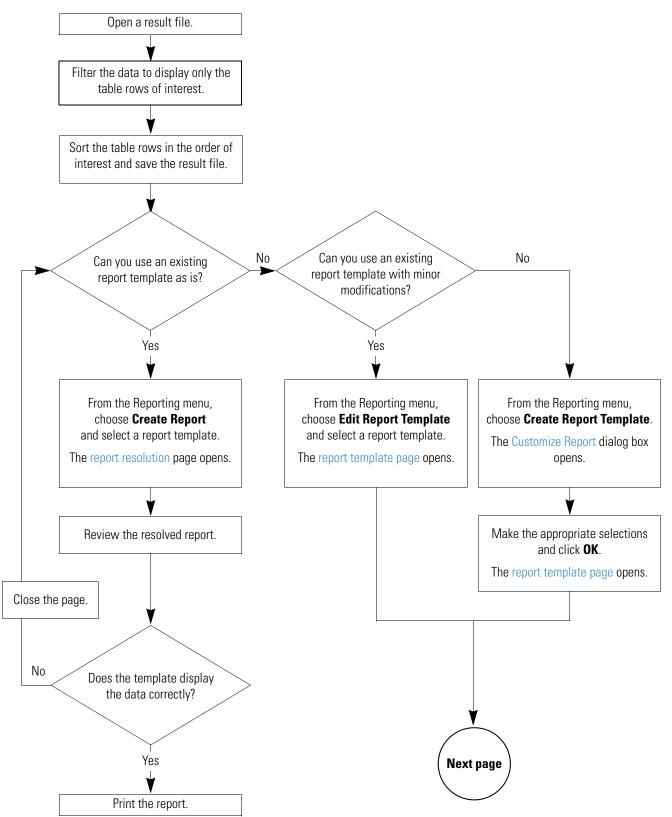
When you open a report template for editing, it opens as a tabbed page with a workspace area on the left

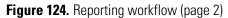
Reporting Workflow

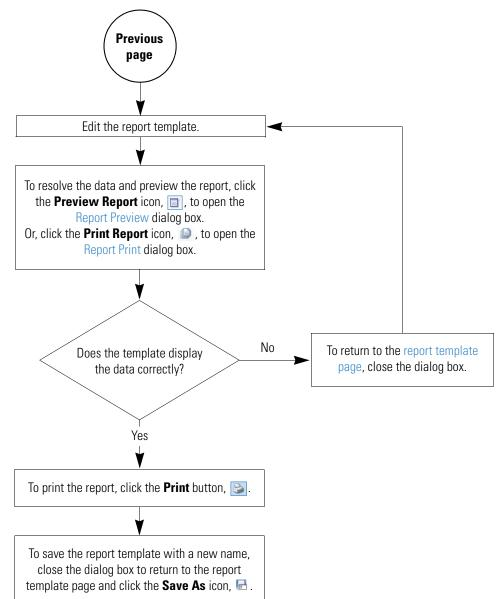
The following flowchart shows the reporting workflow (Figure 123 and Figure 124).

Reporting Workflow

Figure 123. Reporting workflow (page 1)







Generating a Report with an Existing Report Template

You can use one of the report templates provided with the application or one of your own custom report templates to produce reports that display items of interest in a result file.

The application comes with the following eight defined report templates:

- Compounds No Graphs A4.cdReportTemplate
- Compounds with Graphs A4.cdReportTemplate

- Expected Compounds No Graphs.cdReportTemplate
- Expected Compounds per File No Graphs A4.cdReportTemplate
- Expected Compounds per File with Graphs A4.cdReportTemplate
- Expected Compounds with Structures No Graphs.cdReportTemplate
- Expected Compounds with Structures with Graphs A4.cdReportTemplate
- Expected Compounds with GraphsA4.cdReportTemplate

You can find these report templates in the following folder:

C:\Users\Public\Public Documents\Thermo\Compound Discoverer 3.1\Common Templates\ReportTemplates

* To preview and print a report by opening an existing report template

1. Open a result file.

In the application window, the reporting menu commands and the reporting toolbar icons (() become available.

- 2. Determine which main table you want to include in the report and filter the data in this table.
- 3. To select an existing report template, choose **Reporting > Create Report** from the menu bar or click the **Create Report** icon, ().

The Open Report Design Template dialog box opens to the Report Templates folder.

To select an appropriate report template, you must know what data the report template is designed to resolve. Typically, a report template resolves the filtered data from one of the main tables and one or more of the graphs associated with the table. A report template can also resolve data from one or more related tables.

4. Select the appropriate report template and click **Open**.

The report resolution page opens with the thumbnail pane on the right and a report preview on the left.

The tab format for the report resolution page is as follows:

🖲 Report Template Name

As the application resolves the data with the report template, the following icon displays the progress.



When the data is resolved, the progress icon disappears, and the application begins rendering the report pages. The current page/estimated pages box lists the progress.

If the selected template does not contain ReportInfo items, the application displays the pages as it renders them. If the selected template contains a ReportInfo item, the application does not display the rendered pages until it has rendered all of the report pages. ReportInfo items include the time stamp in the upper left and the page number at the center bottom of the defined report templates.

If the report contains too many pages, the application cancels the report generation, and the following message box appears.



- 5. If the application cancels the report generation, repeat step 2 to step 4. But this time, reduce the number of reported items by modifying the filters.
- 6. Review the contents of the report.
- 7. On the report resolution page, click the **Print** icon, 🔄 , in the toolbar to print the report.

The Print dialog box opens.

8. Select the appropriate printer and the page range that you want to print.

The report templates that come with the application default to printing on A4 paper.

- 9. If you are not printing on A4 paper, change the printer setting.
- 10. Click **OK** to print the report.

Related Topics

- Opening, Closing, and Updating Result Files
- Using Result Filters for Data Reduction and Creating Filter Sets

Creating a New Report Template

This topic describes how to create a new report template by using the Customize Report dialog box where you do the following:

- Select the data to be included in the report:
 - Columns of interest in the main table
 - Graphs associated with the main table
 - Columns of interest in any of the related tables
 - Graphs associated with any of the selected related tables
- Change the appearance of the tables:
 - Change the column orientation from left to right to from top to bottom
 - Add Separator lines below the column headings
 - Indent the related tables below the main table
 - Select background colors for the table headers and table rows
- Select the paper type (PaperKind parameter), page orientation, and logo image

See these topics:

- Creating a New Report Template by Using the Customize Report Dialog Box
- Adding, Removing, or Modifying Color Schemes for a Report Template's Tables
- Customize Report Dialog Box Parameters

Creating a New Report Template by Using the Customize Report Dialog Box

To create a report template by using the Customize Report dialog box

- 1. Open a result file.
- 2. Do one of the following:
 - From the menu bar, choose **Reporting > Create Report Template**.
 - In the toolbar, click the **Create a New Report Template** icon, **(**

The Customize Report dialog box opens in front of the New Report Template page.

Oustomize Report - - -Reported Table Compounds -▲ Compounds Columns ġ Structure Name Formula Annotation Source 🗹 Draw Lines 🔲 Transpose Data Indenting Appearance 0.00 Reset Color Scheme Transparent/Transparent Modify -General Settings PaperKind A4 Ŧ Orientation Portrait • Logo Image <u>0</u>K Cancel

Figure 125. Customize Report dialog box

3. In the Reported Table list at the top of the dialog box, select the main table for the report template.

A list of data items for the selected table appears. By default, the Columns list is expanded, and the Graphs and Related Tables lists are collapsed.

- 4. To select the columns for the main table, any of the associated graphs, and any of the main table's related tables, do the following:
 - a. Under the selected table name, click each expand icon, **>**, to open these sections as needed:
 - Columns
 - Graphs
 - Related Tables
 - b. In the expanded sections, select the check box for each column, graph, or related table (and associated columns) that you want to include in the generated reports.
- 5. To add a separator line above the column heading row for each result table, select the **Draw Lines** check box.

(Default layout for the main result table)							
Column 1 label	Column 2 label	Column 3 label	Column 4 label	Column 5 label			
Data text box Data text box Data text box Data text box							
(Related table)							
Column 1 label	Column 2 label	Column 3 label	Column 4 label	Column 5 label			
Data text box	Data text box	Data text box	Data text box	Data text box			

Tip For result tables with a large number of data columns, transpose the columns by selecting the Transpose Data check box. Otherwise, you might need to resize the data columns to fit the page.

- 6. To transpose the tabular data from columns to rows, do the following:
 - a. Select the main result table and the related tables for the report.
 - b. In the data item list, select the table that you want to transpose as follows:
 - To transpose the columns in the main table, select the main table name (the first data item) in the Customize Report dialog box.
 - To transpose the columns in a related table, select the check box to the left of the related table name and click the table name to make sure that it is highlighted in blue.
 - c. Select the Transpose Data check box.

Note Selecting another table clears the Transpose Data check box.

Each selected data column appears as a two-column row in the report template. The first column displays the column heading and the second column displays the data from a table row.

(Default layout for the main table)						
Column 1 label	Column 2 label	Column 3 label				
Data text box	Data text box	Data text box				
(Transposed layout for the main table)						
Column 1 label	Data text box					
Column 2 label	Data text box					
Column 3 label	Data text box					

- 7. To indent a related table, do the following:
 - a. Select the related table in the expanded list of data items.
 - b. In the Indenting box, type the indentation value from **0.00** to **1.00** inch.

(Default layout for the main result table)								
Column 1 label	Column 2 label	Column 3 label	Column 4 label	Column 5 label				
Data text box	Data text box	Data text box	Data text box	Data text box				
				······				
(Related table in	ndented by 1 inch)							
	Column 1 label	Column 2 label	Column 3 label	Column 4 label				
	Data text box	Data text box	Data text box	Data text box				

- 8. To select a color scheme for a report table, do the following:
 - a. In the report item list, select the table of interest, for example, the main table or one of the related tables.
 - b. In the Color Scheme list, select one of the available color schemes.

Each scheme consists of two colors: the first color for the background of the table headers and the second color for the background of the table rows. The default color scheme is Transparent/Transparent.

- 9. Under General Settings, do the following:
 - In the PaperKind list, select the default paper size for the report.
 - In the Orientation list, select the page orientation.
- 10. To apply the settings and close the Customize Report dialog box, click OK.

Your selections appear on the report template page. The tab format for the report template page is as follows:

🔕 Main Result Table

11. To save the template with a different name, click the Save As icon, 🖬 , in the upper left corner of the report template page.

For information about editing the report template, see "Editing an Existing Report Template" on page 450.

When you create a report by using the Customize Report dialog box, the application adds the following additional TextBox design items to the report template: DateTimeInfo in the Page Header section and PageNumberInfo in the Page Footer section.

Figure 126 shows a report template for selected columns in the Compounds per File table and two associated graphs. The default report template uses the Compound Discoverer application icon. You can select a different image and resize its picture container as needed.

Figure 126. Report template with data from the Compounds per File table

- I.	e-and-time stamp to th ected main table name				Logo image's picture container	Default logo image
	CoverPage					
	PageHeader					
	 Compounds per 	File				
	DateTimeInfo					2
Selected	Molecular Weight	RT [min]	Max. # MI	# Adducts	Area	Study File ID
columns	DetailSection_Compound	inds_per_File				
	Molecular Weight	RT [min]	Max. # MI	#Adducts	Area	ĉ
Selected —— graphs	12 10 08 06 04 02 00 00 02 00 □ PageFooter © Reported with Compour □ Appendix			2 0 18 16 14 12 0 00 02 2 2 ageNumberIn	04 06 00 — Adds a page n	
					the report.	

Adding, Removing, or Modifying Color Schemes for a Report Template's Tables

In the Customize Report dialog box, follow this procedure to modify, add, or remove color schemes.

- * To modify the current color scheme or to add or remove color schemes from the list
- 1. In Customize Report dialog box, select the main result table and the related tables for the report.
- 2. Click Modify.

Two color selection lists appear below the Color Scheme list. The list on the left changes the background color for the column headings. The list on the right changes the background color for the data columns.

Color Scheme	Transparent/Transparent	Modify	
	Add	Remove	

3. From the color lists, select one or two background colors, and then click Add.

Oustomize Repo	t	- • ×
Reported Table	Compounds v	
 ✓ Compounds ✓ Columns ✓ Check ✓ Name ✓ Predia 		
Appearance	Draw Lines Transpose Data Indenting 0.00	Reset
Color Scheme	Transparent/Transparent v	Modify
General Settings	Add Remove	
PaperKind	A4 ~]
Orientation	Portrait 🔹]
Logo Image		
	ОК	Cancel

The application displays the effect of the color scheme in the data item list, adds the new color scheme to the Color Scheme list, and activates the Remove button.

② Customize Repo	rt		- • ×
Reported Table	Compounds	-	
 ▲ Columns ✓ Columns ✓ Check ✓ Name ✓ Predia 			۵ ۲
Appearance	Draw Lines Transpose Data Indenting	0.00	Reset
Color Scheme	Turquoise/PaleTurquoise	•	Modify
	Add	Remove	

Note Accepting the settings in the Customize Report dialog box adds the new color scheme to the ColorScheme.xml file that is located in the following folder:

C:\Users\Public\Public Documents\Thermo\Compound Discoverer 3.1\Common Templates\ReportTemplates

If you remove the new color scheme before you click OK at the bottom of the Customize Report dialog box to accept the settings, the application does not add the new color scheme to the ColorScheme.xml file.

- 4. To change the color selection, do one or both of the following:
 - If you do not want to apply the new color scheme to the currently selected table, click **Reset**.

The application undoes the color selections, applies the default color scheme (Transparent/Transparent), and closes the color lists. When you click OK to accept the settings and close the Customize Reports dialog box, the application adds the new color scheme to the ColorScheme.xml file.

• If you do not want to keep the new color scheme, click Remove.

The application undoes the color selections, leaves the color lists open, and removes the color scheme from the Color Scheme list.

Customize Report Dialog Box Parameters

Table 136 describes the parameters in the Customize Report dialog box.

Parameter	Description			
Reported Table	Lists the main tables in the result file.			
Columns	Lists the columns for the selected main table.			
Graphs	Lists the graphs for the selected main table.			
Related Tables	Lists the tables related to the selected main table.			
Columns	Lists the columns for the selected related table.			
Graphs	Lists the graphs for the selected related table.			
Related Tables	Lists the second-level related tables for the selected related table.			
Appearance				
Draw Lines	Specifies whether the application draws a line above the table column headers.			
	Default: Selected			

Table 136. Customize Report dialog box parameters (Sheet 1 of 3)

Parameter	Description				
Transpose Data	Specifies the layout of the data in the result table columns.				
	The default layout (check box cleared) matches the result table layout, with columns displayed from left to right and rows displayed from top to bottom. Select this check box to transpose the columns to rows. Default: Cleared				
	Tip When you select a table item in the data item list, the application automatically clears the Transpose Data check box. For each table that you want to transpose, select the table name and make sure that it is highlighted in blue. Then select the Transpose Data check box.				
Indenting	Specifies the indentation of the selected related table data from the left edge of the page, from 0.00 to 1.00 inch.				
	Default: 0.00 in.				
Color Scheme					
Color Scheme	Specifies the color scheme for the selected table.				
	Each color scheme consists of two colors. The first color is the background of the column headings. The second color is the background of the table rows.				
	You can modify the Color Scheme list by adding or removing color schemes.				
	Default: Transparent/Transparent				
	Note Accepting the settings in the Customize Report dialog box adds the new color schemes to the ColorScheme.xml file that is stored in the same folder as the common report templates.				
General Settings					
PaperKind	Specifies the size of the paper for printing the report. Select the appropriate paper size before sending the report to the printer.				
	Default: A4				
Orientation	Specifies the orientation of the report, either Portrait or Landscape.				
	Default: Portrait				

Table 136. Customize Report dialog box parameters (Sheet 2 of 3)

Description
Specifies the logo image to appear by default in the upper right corner of each report page.
The default size of the picture container for the logo is 1.823×0.492 in. (width × height). When the selected image is larger than the picture container, the container clips the image. You can edit the properties of the picture container in the report template.
Resets the color scheme to the default scheme.
Opens two color selection lists.
Selecting colors in one or both of the color selection lists below the Color Scheme list activates this button.
Applies the new color scheme to the selected table and adds the new color scheme to the Color Scheme list.
Removes the selected color scheme from the Color Scheme list.
Applies the selected settings to the new report template.
Cancels your selections and closes the dialog box.

Table 136. Customize Report dialog box parameters (Sheet 3 of 3)

Editing an Existing Report Template

Use the report template page to modify a report template. The report template page shows the design items that you selected using the Customize Report dialog box or the items in the existing report template that you selected. Some of the design items appear as containers (a rectangular box) where you can add text, images, or data graphs.

Note You can open more than one report template page in the application window.

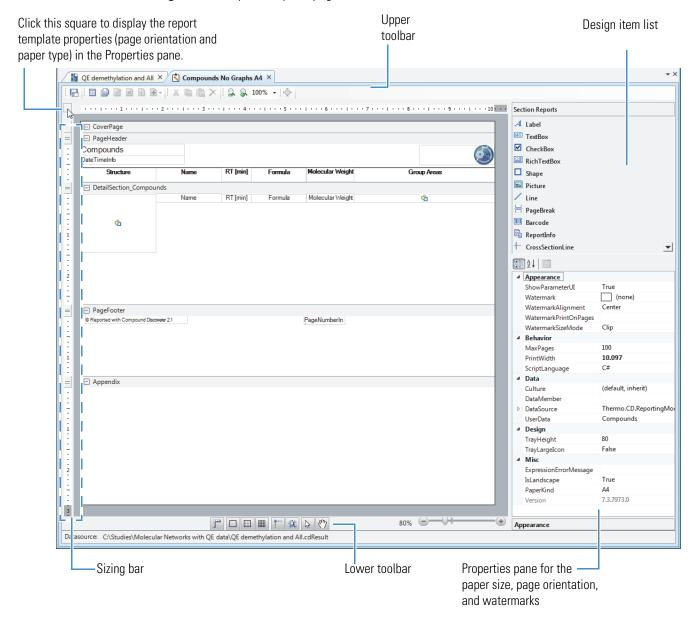
To modify a report template, see these topics as needed:

- Opening a Report Template
- Changing a Report Template's Logo Image
- Changing a Report Template's Time Stamp
- Adding Design Items to a Report Template
- Adding Main Table Columns to a Report Template
- Adding Data Graphs to a Report Template

- Adding Related Table Columns to a Report Template
- Editing the Properties of Subreport Columns in a Report Template
- Moving a Subreport Column to a Report Template's Main Table
- Adding Page Breaks to a Report Template
- Deleting a Pair of Workspace Sections on the Report Template Page
- Resizing a Report Template's Sections
- Adding Design Items to a Report Template
- Adding Main Table Columns to a Report Template
- Adding Data Graphs to a Report Template
- Adding Related Table Columns to a Report Template
- Editing the Properties of Subreport Columns in a Report Template
- Moving a Subreport Column to a Report Template's Main Table
- Modifying the Properties of a Report Template Design Item
- Adding Page Breaks to a Report Template

Figure 127 shows the report template page.

Figure 127. Report template page



Related Topics

- Editing Tools for the Report Template Page
- Selecting the Paper Type, Print Width, Page Orientation, and Watermark for a Report Template

Opening a Report Template

To open a report template

- 1. Open a result file.
- 2. From the application window, do one of the following:
 - a. Choose **Reporting > Create Report Template** from the menu bar, or click the **Create Report** icon, (**b**), in the toolbar.

The Customize Report dialog box opens.

b. Make the appropriate selections and click **OK**.

The report template page opens as a tabbed document. The tab format is as follows:

🔕 Main Result Table Name

-or-

a. Choose **Reporting > Edit Report Template** from the menu bar, or click the **Edit Report Template** icon, (1), in the toolbar.

The Open Report Design Template dialog box opens.

b. Select a report template and click **Open**.

The report template page opens as a tabbed document. The tab format is as follows:

S Existing Report Template Name

Changing a Report Template's Logo Image

To change the logo image in a report template

1. Open the report template.

The report template opens as a tabbed page in the application window.

2. Select the logo image container.

The picture properties appear in the Properties pane.

Note The Compound Discoverer icon is the default logo for the common templates and the templates that you create with the Customize Reports dialog box.

3. In the Data area, click the browse icon,, to the right of the Image property. You might have to click the row to make the browse icon appear.

Image System.Drawing.Bitmap

....

The Open dialog box opens with a setting of All image files for the file type.

4. Browse to the folder where you stored the logo of interest, select the logo, and click **Open**.

The selected image appears in the container.

5. Modify the Layout properties as appropriate.

Changing a Report Template's Time Stamp

* To change the format of the date-and-time stamp

1. Open the report template.

The report template opens as a tabbed page in the application window.

2. Select the DateTimeInfo item.

The properties for this design item appear in the Properties pane.

3. Under Appearance, click the **OutputFormat** box to make the browse icon appear. Then click the browse icon.

The OutputFormat dialog box opens (Figure 128).

Figure 128. OutputFormat dialog box

<u>C</u> ategory: General Number	Sample 05-Apr-1998 2:30	
Currency Date Time	dd-MMM-yyyy h:mm	
Percentage Custom	MMMM-yy MMMM-yyy MMMM d, yyyy M/d/yyy h:mm tt M/d/yyy h:mm tt M/d/yyy h:mm M/d/yyy h:mm M/d/yyy h:mm dd-MMM-yyyy h:mm dd-MMM-yyyy h:mm:ss	

- 4. Select the format of interest.
- 5. Click **OK** to accept the setting.

Adding Design Items to a Report Template

You can add any of the design items in the Section Reports pane to a report template.

To add a design item to the template

1. Open the report template.

The report template opens as a tabbed page in the application window.

2. From the Section Reports pane to the right of the workspace area, drag a design item to the appropriate location on a workspace section of the page.

Related Topics

• Section Report Items for a Report Template

Adding Main Table Columns to a Report Template

Use the Add Field shortcut menu command or the Add Items icon in the upper toolbar to add data fields from a result table to the report template.

To add a main table column to the report template

1. Open the report template.

The report template opens as a tabbed page in the application window.

- 2. Depending on the orientation of the table columns, do the following:
 - If the table columns are arranged horizontally from left to right, go to step 3.
 - If the table columns are arranged vertically from top to bottom, go to step 5.
- 3. To add a main table column to a column set that is arranged horizontally, do either of the following:
 - To place the new column to the right of the current column set, select the **PageHeader** bar.
 - To place the new column to the right of a specific column, select the column heading.
- 4. Do one of the following:
 - Right-click the PageHeader bar or a specific column heading, and then choose Add Field > *Data Column* from the shortcut menu.

-or-

- a. Click the PageHeader bar or a specific column heading.
- b. In the toolbar, click the **Add Items** icon, **e**, to open a list of data column selections.

The available items in the list include the unused data columns in the current main table. The data columns that are already in the template are unavailable and grayed out.

c. Select an available data column from the list.

The new data column appears to the right of the selected column or the current column set. If there is a gap to the right of the selected column or to the right of the column set, the new column fills the gap. If there is no space to the right of the selected column or to the right of the column set, the new column shares the space with the selected column or the last column in the column set.

- 5. To add a main table column to a column set that is arranged vertically, do the following:
 - a. Use the sizing bar to display all of the data column rows.
 - b. Right-click the Label column (heading) of the two-column row that is above where you want to add the new two-column data row, and choose Add Field > Data Column.

The new two-column data row appears below the selected two-column row.

Adding Data Graphs to a Report Template

* To add a data graph that is associated with the main table to the template

1. Open the report template.

The report template opens as a tabbed page in the application window.

- 2. Do one of the following:
 - Right-click the DetailSection_Main_Table_Name bar and choose the Add Field > Data Graph of interest from the shortcut menu.

-or-

- a. Click the **DetailSection_***Main_Table_Name* bar.
- b. In the toolbar, click the **Add Items** icon, **e**, to open a list of data graphs and related table selections.

The available items in the list include the data graphs associated with the current main table and the related tables for the current main table.

c. Select an available data graph from the list.

When the table columns are positioned from left to right, the data graph appears below the table columns and to the right of any existing data graphs.

Adding Related Table Columns to a Report Template

To add a column from a related table to a report template

1. Open the report template.

The report template opens as a tabbed page in the application window.

2. Do one of the following:

CoverPage							
PageHeader							
Expected Com	npounds per Fi	le					
DateTimeInfo							
Parent Compound	Formula	Мо	lecular V	Veight	Compositi Change		RT [min]
DetailSection PageFooter @ Reported with C Appendix	Add Field Align Fields Transpose Data Field Insert Section Reorder Group Paste		52 Spectr romatog pected C put Files	rum gram Trace		Checked	
	Delete Properties		· · ·	lated Str			Ion Charge
							Molecular Weight m/z ΔMass [Da] ΔMass [ppm] RT [min]

 Right-click the DetailSection_Main_Table_Name bar and choose Add Field > Related Table Name > Column of Interest.

-or-

- a. Select the **DetailSection_***Main_Table_Name* bar.
- b. In the toolbar, click the **Add Items** icon, **I**, to open a list of data graphs and related table selections.

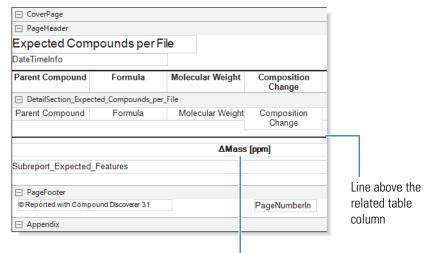
The available items in the list include the data graphs associated with the current main table and the related tables for the current main table.

- c. Select an available related table from the list, and then select a table column.
- 3. If the DetailSection is collapsed, click the expand icon to open the section.

DetailSection_Expected_Compounds_per_File
 Expand icon

The added table column appears at the bottom of the DetailSection (data area). The application automatically adds a line above the column heading (Figure 129).

Figure 129. Expected Compounds per File table with the Δ Mass [ppm] column from the related Expected Features table



Column from the related table

Editing the Properties of Subreport Columns in a Report Template

To edit the properties of subreport columns

1. Open the report template.

The report template opens as a tabbed page in the application window.

- 2. To open the subreport editor area, do one of the following:
 - Double-click the **Subreport_***Related_Table_Name* box.

-or-

a. Select the **Subreport_***Related_Table_Name* box.

The Edit Sub-Report icon, 📝, becomes available.

b. Click the **Edit Sub-Report** icon.

The TextBox design item for the related table column appears in a separate section. The item's container is sized to the full page width (Figure 130).

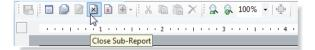
CoverPage				1
PageHeader				
Expected Comp	ounds per Fil	e		
DateTimeInfo				
Parent Compound	Formula	Molecular Weight	Composition Change	
DetailSection_Expect	ed_Compounds_per_l	File	-	
Parent Compound	Formula	Molecular Weight	Composition Change	
	ΔΜ	ass (ppm)		
Subreport_Expected_f	Features			
PageFooter				
© Reported with Compo	und Discoverer 3.1		PageNumberIn	
Appendix			···· 🕸 📐 🖑	
	1 2 1		. 4	
🖃 🚺 DetailSection_Exp	ected_Features			
			ΔMass [ppm]	
olumn heading for Mass [ppm] colum			Subreport section	with the containe
			the Δ Mass [ppm] full page width	data item sized to

Figure 130. Report template page with the subreport section open

3. To change the properties of the TextBox design item in the subreport section, select it.

The properties for the selected design item appear in the properties pane to the right of the workspace.

- 4. Make changes as necessary in the properties pane, or click the property dialog link below the properties pane to open the TextBox dialog box and make similar changes.
- 5. To close the subreport section, click the **Close Sub-Report** icon in the report designer toolbar.



Moving a Subreport Column to a Report Template's Main Table

To move a subreport column up to the set of main table columns *

1. Open the report template.

The report template opens as a tabbed page in the application window.

2. Resize the column heading container for the subreport column. Then, move it to an appropriate location in the set of main table columns.

- 3. Resize the Subreport_*Related_Table_Name* container. Then, move it to the appropriate location in the Detail Section.
- 4. In the related table's Detail Section, resize the data item's container.

Figure 131 shows a report template for the Expected Compounds per File table with an additional column from the related Expected Features table.

Figure 131. Related table column moved up to the main table

				Сс	olumn h	eading for th	ne
				re	lated ta	ble column	
CoverPage							
🖃 PageHeader							
Expected Co	ompounds	per File	Э				
DateTimeInfo							
Parent Fo Compound	ormula		lecular Neight	ΔMa	ss (ppm)	Composition Change	
DetailSection_E	xpected_Compou	unds_per_F	ile D.	Th	00	_	
Parent Fo Compound	rmula	Molecular		Subrep		Composition	
			0				
PageFooter							
© Reported with Co	ompound Discover	er 3.1			P	ageNumberIn	
Appendix							
		ł				🛱 🏷 🖑	
(1,1,1,1,1,1,1,1,1,1,1,1,1,1,1,1,1,1,1,		2	· · · 3		4		
DetailSection	- Expected Feat	uror.			_		
ΔMass [ppm]	n_Expected_Featu	1103					
d							
				c	ubrono	t containar	moved to the
						le's Detail S	
				11		ies Delali S	ection
Detail Sec	ction for the	related	l table	with	а		
	ontainer for				-		

Modifying the Properties of a Report Template Design Item

On the report template page, you can edit the position and size of a design item by using the mouse or the properties pane on the right. You can also edit other properties for a design item from the properties pane or the specific dialog box for the design item.

To edit a design item

1. Open the report template.

The report template opens as a tabbed page in the application window.

2. Select the design item.

The properties for the selected design item appear in the properties pane, at the bottom right of the report template page. For information about the properties of each design item, see "Property Settings for the Sections and Items in a Report Template" on page 474.

- 3. Do any of the following:
 - To move the design item to another position, do either of the following:
 - Move the item by dragging it (or you can use the arrow keys on the keyboard).
 - In the Properties pane, under Layout, expand Location, and then change the X and Y values.
 - To resize the container for the design item, do either of the following:
 - Resize the item by dragging the handle points of the container.
 - In the Properties pane, under Layout, expand Size, and then change the Width and Height values.
 - To change other properties for the design item, do either of the following:
 - Modify the properties in the Properties pane.
 - Click the **Property Dialog** link at the bottom right of the report template page, and then modify the property in the item-specific dialog box.

Note For the RichTextBox design item, in addition to the Property Dialog link, you can click the Load File link to load text from a file (see "Populating a Rich Text Box by Loading a File" on page 487).

Adding Page Breaks to a Report Template

- * To add a page break between reported result table rows
- 1. Open the report template.

The report template opens as a tabbed page in the application window.

2. On the report template page, drag the PageBreak design item from the Section Reports pane to the bottom of the Detail Section of the report template.

Deleting a Pair of Workspace Sections on the Report Template Page

Note The report template page pairs these sections together:

- Cover Page and Appendix
- Page Header and Page Footer

When you select one of the paired sections to delete, the application removes both sections. You cannot delete one section without deleting the other, and you cannot delete the Detail Section section.

To delete a pair of workspace sections

1. Open the report template.

The report template opens as a tabbed page in the application window.

- 2. Select one of the workspace sections by clicking the section header or by clicking within the section area.
- 3. From the shortcut menu, choose **Delete**.

Resizing a Report Template's Sections

Use the sizing bar to the left of the workspace to resize each workspace section vertically, except for the Appendix section.

Tip Make sure to enlarge the workspace section enough to hold all of the items that you want to add to that section of the template.

The size of a section on the report template page is not necessarily the same as its size in the generated report.

✤ To vertically resize a workspace section

1. Open the report template.

The report template page opens as a tabbed page in the application window.

- 2. Do any of the following:
 - For the CoverPage, PageHeader, DetailSection, and PageFooter sections, do the following
 - In the sizing bar, drag the sizing handle.
 - To vertically enlarge a workspace section, drag down the handle that is aligned with the header of the subsequent section. To reduce a workspace section, drag the handle up.
 - To enlarge the Appendix section, drag it down by the bottom edge of the report template page. To reduce this workspace section, drag the bottom edge up.

Related Topics

• Workspace Sections and Sizing Bar on the Report Template Page

Editing Tools for the Report Template Page

For information about the editing tools for the report template page, see these topics:

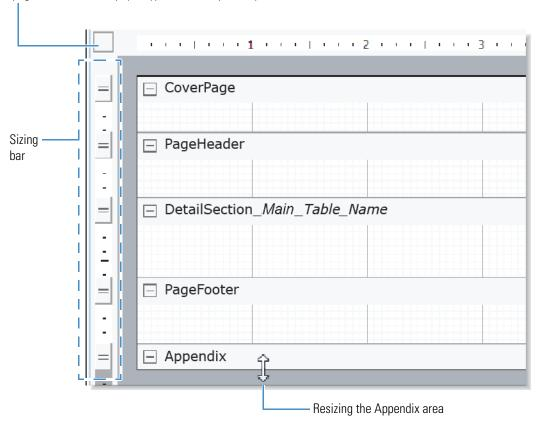
- Workspace Sections and Sizing Bar on the Report Template Page
- Report Template Page Toolbars
- Report Template Page Shortcut Menu
- Section Report Items for a Report Template
- Property Settings for the Sections and Items in a Report Template
- Opening the Property Dialog Boxes for the Report Template's Design Items
- Populating a Rich Text Box by Loading a File

Workspace Sections and Sizing Bar on the Report Template Page

Figure 132 shows the five workspace sections, the sizing bar on the left, and the icon that opens the page orientation and paper type options in the upper left corner.

Figure 132. Workspace sections, sizing bar, and icon that opens the page size options

Click this square to display the report template properties (page orientation and paper type) in the Properties pane.



By default, the workspace on the report template page has five sections. Table 137 lists these sections, from top to bottom.

Workspace section	Description					
Cover Page	Appears as the first page of a report with the following property selections:					
	• True—For the Visible parameter in the Behavior section of the properties pane.					
	Default: False					
	• After—For the NewPage parameter in the Data section of the properties pane.					
	Default: After (See "NewPage" on page 482.)					
	To display the Cover Page section without a page break between it and the next section, select True for the Visible parameter and None for the New Page parameter.					
	Use this section to add nonrepeating information, such as the report title, date-and-time stamp, and company logo.					
Page Header	Adds design items to the top of each report page. The standard templates include a Label design item with the main table name, a TextBox design item with a time stamp, and a Picture design item with a company logo. The column headings (Label design items) appear here when you add table columns to the report.					
Detail Section (concatenated with the selected <i>Main_Table</i> name)	Adds data from the result file, such as the repeating items (TextBox design item) of a main table or the data graphs for each table row.					
Page Footer	Adds information to the footer of each report page, for example, the page number.					
	Report templates created with the Customize Reports dialog box automatically include a page number at the bottom of each page. The page number is a TextBox design item.					

Table 137. Default workspace sections (Sheet 1 of 2)

Table 137. Default workspace sections (Sheet 2 of 2)				
Workspace section	Description			
Appendix	Adds information to an appendix section of the report.			
	An Appendix section appears after the last page of a report, with the following property settings:			
	• True—For the Visible parameter in the Behavior section of the properties pane			
	Default: False			
	• Before—For the NewPage parameter in the Data section of the properties pane			
	Default: Before			
	To display the Appendix section without a page break between it and the previous section, select True for the Visible parameter and None for the NewPage parameter.			

Table 137. Default workspace sections (Sheet 2 of 2)

Report Template Page Toolbars

Table 138 describes the toolbars on the report template page.

For more information about the report template page, see "Editing an Existing Report Template" on page 450.

Table 138. Toolbars on the report template page (Sheet 1 of 4)

lcon	Description
Top toolbar	
	Save Active Item—Saves the report template using the same file name.
	By default, the template file name is the same name as the main table that you selected in the Customize Report dialog box (see "Creating a New Report Template" on page 442).
Ξ	Save As—Saves the report template using a different file name.
	Preview—Opens the Report Preview dialog box (see "Previewing and Printing a Report" on page 489).
	Print—Opens the Report Print dialog box.

con	Description					
2	Edit Sub-Report—Enlarges a selected subreport so that you can edit it.					
	Selecting a subreport item activates this icon. When you click this icon, the report designer opens the subreport in a separate Detail Section workspace section. You can zoom in on this temporary section or zoom out of it. Increasing the size of this temporary section does not affect the report template page.					
	Note Related tables that you select in the Customize Report dialog box appear as subreports on the report template page.					
×	Close Sub-Report—Closes the separate subreport workspace section.					
	Clicking anywhere in the separate subreport workspace activates this icon.					
	Align Columns—Aligns the Label (column heading) and Textbox (data) containers for the selected column or columns.					
	Selecting a report column activates this icon.					
.	Add Items—Opens a list of items that you can add to the currently selected section of the report template.					
	Clicking within a workspace section or the section header activates this icon. The list of items varies depending on the selected section:					
	• CoverPage and DetailSection sections: You can add related table columns or graphs that are not currently in the template					
	• PageHeader section: You can add main table columns that are not currently in the template. The column heading appears ir the PageHeader section as a Label design item, and the container for the column data appears in the DetailSection as a TextBox design item.					
x	Cut—Deletes the selected item without confirmation.					
Ē	Copy—Copies the selected item.					
a	Paste—Pastes the selected item.					

Table 138. Toolbars on the report template page (Sheet 2 of 4)

lcon	Description				
2	Zoom Out—Reduces the magnification of the page.				
\$	Zoom In—Increases the magnification of the page.				
100% -	Magnification box—Displays the magnification percentage.				
١	Actual Size—Displays the page at 100% magnification.				
	Changing the magnification by using the Zoom In and Zoom Out icons or by typing a value in the magnification box activates this icon.				
Bottom toolbars					
	Dimension Lines—Displays dimension lines (—1.0 in—) as you resize a design item by using the mouse.				
	Hide Grid—Clears the grid on the page.				
	Show Dots—Shows the main grid lines and the small dots within the grid.				
	Show Lines—Shows the main grid and the smaller lines within the grid.				
** ***	Snap Lines—When you move a design item on the page, blue alignment lines appear. When the selected design item (Item 1 below) is horizontally aligned with another design item, two vertical lines bracket the aligned items. When the selected design item is vertically aligned with another design item, two horizontal lines bracket the aligned items.				
	Item 1 Item 3				
	Snap to Grid—When you move an item on the page, this mode automatically snaps it to the smaller grid lines.				
×	Select Mode—Use this mode to select items on the page.				
	To select multiple items, press the SHIFT key while you select the items.				

Table 138. Toolbars on the report template page (Sheet 3 of 4)

lcon	Description
٣ ٦	Pan Mode—When the page is zoomed in, use this mode to move to a different part of the page.
100%	Magnification bar—Move the slider to the left to zoom out and to the right to zoom in. The magnification percentage appears to the left of the slider.

Table 138. Toolbars on the report template page (Sheet 4 of 4)

Report Template Page Shortcut Menu

Follow these procedures as needed.

* To open the shortcut menu for the report template page

Right-click the report template page.

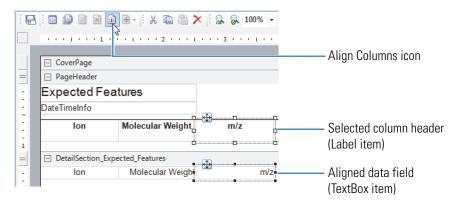
* To add a design item to a section of the report

Right-click the section bar within a section area and choose Add Field > Item of Interest.

✤ To align a column heading to its associated data field

1. In the PageHeader section, right-click the column heading (Label design item) that you want to align with its associated data field (TextBox design item).

The Align Columns icon becomes available.



2. From the shortcut menu, choose Align Fields.

Tip You can also select the column heading, and then click the **Align Columns** icon, i , in the report template page toolbar.

- To transpose the data fields from columns to rows or from rows to columns
- 1. Right-click a column heading (Label) or a column data field (TextBox) and choose **Transpose Data Fields**.

2. To undo the change, if necessary, right-click a column heading (Label) or a column data field (TextBox) and choose **Transpose Data Fields**.

✤ To add a border to a design item

1. Right-click the design item in the workspace and choose Format Border.

The Format Border dialog box opens (Figure 133).

Figure 133. Format Border dialog box with the selection of a coral, double line border

Format Border - ActiveReports					
Presets	Line Styles	Preview Click on diagram below or use presets to edit borders Image: Click on the set of			

2. Click one of the icons in the Presets area or select the line style in the Line Styles area, and click the appropriate sides of the square in the Preview area to set up the border.

Table 139 describes the shortcut menu commands for the report template page.

Table	139.	Report	template	page	shortcut	menu	(Sheet 1	of 2)

Command	Description			
Add Field	Adds the item that you choose to the selected workspace section.			
Align Fields	Aligns a column header or subreport with the associated data.			
	This command becomes available when you select a column header or a subreport.			
Transpose Data Fields	Transposes the data from columns to rows or from rows to columns.			
	This command becomes available when you select a column header.			
	on command is available if the template does not already include a section or a Page Header/Footer section.			
Insert Section > Report Header/Footer	Inserts the ReportHeader and ReportFooter sections.			
Insert Section > Page Header/Footer	Inserts the PageHeader and PageFooter sections.			

Command	Description
Сору	Copies the selected item.
Paste	Pastes the selected item.
Cut	Removes the selected item without confirmation.
Delete	Deletes the selected item after you click OK in the confirmation dialog box.
Bring to Front	Moves the selected item to the front, on top of other surrounding items.
Send to Back	Moves the selected item to the back, beneath all other surrounding items.
Format Border	Opens the Format Border dialog box where you can change an item's border layout, line style, and color.
Properties	Highlights the (Name) property in the properties pane of the report template page.

Table 139. Report template page shortcut menu (Sheet 2 of 2)

Section Report Items for a Report Template

The Section Reports pane to the right of the workspace on the report template page contains all of the different design items that you can add to the report template.

To add a design item to a report template

1. Drag the item of interest to the appropriate location on the report template page.

Some of the items appear as containers (boxes) where you can add text or images.

- 2. Select the design item to open its properties below the Section Reports pane.
- 3. Edit the item's properties as necessary.

Table 140 describes the design items in the Section Reports pane.

Table 140. Design items in the Section Reports pane (Sheet 1 of 3)

Design item	Description
Label	A text label that is usually used as a header or title. The column headers are Label design items.
TextBox	A text box that is usually used to group multiple items together. The application uses the TextBox design item to display the repeating tabular data in the result file.
	IMPORTANT Use the Rich Text Box item rather than the TextBox item to add text to the report's cover page.

Design item	Description
CheckBox	A check box that you can select or clear.
RichTextBox	A text box that you can populate by typing text in the box or by loading text from a file. Use the Load File link below the properties pane on the right to select and open a text file.
Shape	A geometric shape such as a rectangular or square box (with either square or rounded corners), an ellipse, or a circle. Style Rectangle Rectangle Ellipse RoundRect RoundRect
	Tip When you add this item, by default, it appears as a rectangular box with square corners. To change to a different shape, modify the Style property under Appearance in the properties pane.
Picture	A container for a graphic.
Line	A straight line.
PageBreak	A break to push the subsequent content to the next page. To place reported table rows on separate pages, add a page break at the bottom of the DetailSection.
Barcode	A bar code. For information about setting up the Barcode properties, refer to the <i>ActiveReports User Guide</i> on the company website for Grape City.

Table 140. Design items in the Section Reports pane (Sheet 2 of 3)

Design item	Description
ReportInfo	A variable that the application automatically replaces with real-time data in the generated report. Use the ReportInfo design item to show the current page number or to add a date-and-time stamp. Select the page number or date-and-time stamp from the FormatString list under Appearance in the Properties pane.
	FormatString Page {PageNumber} of {PageCount} on {RunDateTime} Page {PageNumber} of {PageCount} {RunDateTime:} {RunDateTime:M/d} {RunDateTime:M/d} {RunDateTime:M/d/yyy} {RunDateTime:M/d/yyy} {RunDateTime:M/d/yyy} {RunDateTime:M/d/yyy} {RunDateTime:MM/dd/yy} {RunDateTime:MM/dd/yy} {RunDateTime:d-MMM} {RunDateTime:d-MMM-yy} {RunDateTime:MMMM-yy} {RunDateTime:MMMM-yy} {RunDateTime:MMMM-yy} {RunDateTime:MMMM-yyy} {RunDateTime:MMMM-yyy} {RunDateTime:MMMMM-yyy} {RunDateTime:MMMM d, yyyy} {RunDateTime:M/d/yy h:mm tt} {RunDateTime:M/d/yy h:mm} {RunDateTime:M/d/yy h:mm}
CrossSectionLine	A line that can span across multiple workspace sections on the report template page.
	Note You cannot add this item to the Detail Section workspace section. However, you can add it to another section (for example, the Page Header section) and have it span across the Detail Section section.
CrossSectionBox	A box that can span across multiple workspace sections on the report template page.
	Note You cannot add this item to the Detail Section workspace section. However, you can add it to another section (for example, the Page Header section) and have it span across the Detail Section section.

Table 140.Design items in the Section Reports pane (Sheet 3 of 3)

Property Settings for the Sections and Items in a Report Template

On the report template page, the properties pane to the bottom right of the workspace includes all of the property settings that you can use to format a workspace section or an object (design item) in the report template. The settings vary depending on the selected workspace section or design item.

If it is available, click the expand icon, \triangleright , to open the settings for a particular property or the collapse icon, \blacktriangle , to close the settings.

The properties pane contains these property groups, from top to bottom:

- Appearance Properties
- Behavior Properties
- Data Properties
- Design Properties
- Layout Properties
- Miscellaneous Properties
- Opening the Property Dialog Boxes for the Report Template's Design Items
- Populating a Rich Text Box by Loading a File

IMPORTANT The Summary properties are not functional.

Appearance Properties

Table 141 describes the Appearance properties.

Property	Description
AnchorBottom	(For the Line design item only) Specifies whether the line is anchored to the bottom of the workspace section.
	 Selections: False—Does not anchor the line to the bottom of the workspace section. True—Anchors the line to the bottom of the workspace section.
Alignment	(For the Label, TextBox, ReportInfo, and Barcode design items) Specifies the horizontal alignment of the text within the container
	(For the Barcode design item) Specifies the horizontal alignment of the caption text that is associated with the bar code. You enable the caption text by setting the CaptionGrouping and CaptionPosition properties.
BackColor	Specifies the background or fill color.
BarHeight	(For the Barcode design item only) Specifies the height of the bar code.
CaptionGrouping	(For the Barcode design item only)
	 Selections: False—Does not enable a text caption to be associated with the bar code. True—Enables a text caption.
CaptionPosition	(For the Barcode design item only) Sets the position of the caption container relative to the bar code symbol.
	B25
	 Selections: None—Hides the caption. Above—Sets the position above the symbol. Below—Sets the position below the symbol, as in the figure above.
CharacterSpacing	(For the Label and TextBox design items) Specifies the spacing between the characters in the text, in points.
	Default: 0

Table 141. Appearance properties in the Properties pane (Sheet 1 of 3)

Property	Description
ClassName	Specifies the name of the class for a particular format.
	Default: Normal
Font	(For the Label, TextBox, and CheckBox design items) Specifies the name of the font and other font characteristics such as the style, size, effects, and script. Clicking the browse icon,, opens the Font dialog box where you can specify the font characteristics. Clicking the expand icon, ▷, expands the settings.
Font >	Specifies the GDI character set to use. For a list of valid values,
GdiCharSet	refer to the GdiCharSet Property in the Microsoft Developer Network (MSDN) Library.
Font >	Specifies that the font is derived from a GDI vertical font.
GdiVerticalFont	
ForeColor	(For the Label, TextBox, and CheckBox design items) Specifies the font color.
FormatString	(For the ReportInfo design item only) Specifies the format of the generated content as a page number or a date-and-time string.
LineColor	Specifies the color of a line or border.
LineSpacing	Specifies the spacing between multiple lines of content, in points.
LineStyle	Specifies the style of a line or border.
LineWeight	Specifies the thickness of a line or border, in pixels.
NarrowBarWidth	(For the Barcode design item only) Specifies the width of the narrow bars in the bar code (a value of 1.0 equals 0.864 points).
	Tip At a thicker width for the narrow bars, the entire bar code might be too large for the container. In this case, enlarge the size of the container to see the entire bar code.
NWRatio	(For the Barcode design item only) Specifies the ratio of the width of the wide bars relative to the width of the narrow bars in the bar code. The larger the ratio, the thicker the wide bars appear.
	Tip At a thicker width for the wide bars, the entire bar code might be too large for the container. In this case, enlarge the size of the container to see the entire bar code.

Table 141. Appearance properties in the Properties pane (Sheet 2 of 3)

Property	Description
OutputFormat	(For the TextBox design item only) Specifies the format settings for custom content, or for a number, currency, date, time, or percentage. Clicking the browse icon,, for this property opens the OutputFormatDialog box (see Figure 128 on page 454) where you can change the settings.
	Do not change the OutputFormat settings for data fields from a result table column.
PictureAlignment	(For the Picture design item only) Specifies the alignment of the selected image with respect to the container. For proper alignment, the container must be larger than the image.
	Selections: TopLeft, TopRight, Center, BottomLeft, and BottomRight
QuietZone	(For the Barcode design item only) Specifies the left, right, top, and bottom margins of the quiet zone for the bar code, in inches.
Rotation	(For the Barcode design item only) Specifies the rotation of the bar code within the container.
	Tip At a rotation of 90 or 270 degrees, the entire bar code might be too large for the container. In this case, enlarge the size of the container to see the entire bar code.
Style	Specifies the format settings such as color, alignment, font, geometric shape, or bar code properties.
SupplementOptions	(For the Barcode design item only) Specifies the supplement options (2- or 5-digit add-ons for EAN/UPC bar codes).
TextJustify	(For the Label and TextBox design items) Specifies how to distribute the text when you set the Alignment property to Justify
VerticalAlignment	(For the Label and TextBox design items) Specifies the vertical alignment of the text within the container.
	Selections: Top, Middle, and Bottom
VerticalText	 Specifies the a vertical alignment for the text. False—Does not render the text according to the vertical layout settings. True—Renders the text according to the vertical layout settings.
WrapMode	Specifies that a long line of text wraps to the beginning of the next line to fit in the container.

Table 141. Appearance properties in the Properties pane (Sheet 3 of 3)

Behavior Properties

Table 142 describes the Behavior properties.

Table 142. Behavior properties in the Properties pane (Sheet 1 of 3)

Behavior property	Description
Angle	Specifies the slope of the text within the container, in degrees.
AutoReplaceFields	 (For the RichTextBox design item only) Specifies whether the data in the container is automatically replaced with the data from the data source as specified by the Data Field property selection. False—Does not automatically replace the fields of the object with the fields in the data source that are assigned to the current workspace section. True—Automatically replaces the fields of the object with the fields in the data source that are assigned to the fields in the data source that are assigned to the current workspace section.
AutoSize	(For the Barcode design item only) Specifies whether the barcode design item stretches to fill its container. Selections: True or False
CanGrow	Specifies whether the container or section can increase in height to fit its contents.
	Selections: True or False
CanShrink	Specifies whether the container or section can decrease in height to fit its contents.
	Selections: True or False
CheckAlignment	(For the CheckBox design item only) Specifies the alignment of the check box in the container.
Checked	(For the CheckBox design item only) Specifies whether the CheckBox design item appears with or without a check mark in the report.
	Selections:False—Shows the check box selected.True—Shows the check box cleared.
CheckSumEnabled	(For the Barcode design item only) Specifies whether the application computes and includes a checksum in the bar code.
	Selections: True or False

Behavior property Description		
ColumnDirection	(For the Detail Section workspace section only) Specifies whether to display the data columns in the down-and-across direction or the across-and-down direction, for a multi-column (newspaper-style) report.	
	Selections: DownAcross or AcrossDown	
Enabled	(For the PageBreak design item only) Specifies whether the PageBreak design item is enabled.	
	Selections: True or False	
KeepTogether	(For the Detail Section and Appendix workspace sections only) Specifies whether the contents of the current section prints on a single page. This property does not shrink items to fit; rather, it acts like a page break between reported table rows.	
	Selections: True or False	
MultiLine	(For the Label, TextBox, ReportInfo, and RichTextBox design items only) Specifies whether the report template displays only the content that fits on one line or displays all the lines that fit in the container.	
	Selections: True or False	
PrintAtBottom	(For the Appendix workspace section only) Specifies where the design items in the Appendix section are printed.	
	 Selections: False—Places the Appendix design items immediately after the DetailSection and before the page footer information. True—Places the Appendix design items at the bottom of the current page just above the page footer information. 	
RepeatToFill	(For the Detail Section workspace section only)	
	False—Does not repeat content to fill the report page.	
	True—Repeats content to fill the report page.	
RightToLeft	((For the Label, TextBox, and ReportInfo design items) Specifies whether the text is aligned with the right side of the container. Also supports locales that use right-to-left fonts.	
	Selections: True or False	

Table 142. Behavior properties in the Properties pane (Sheet 2 of 3)

Behavior property	Description		
ShrinkToFit	Specifies whether the font size of the text within the selected container shrinks to fit the container. If the WrapMode property under Appearance is set to WordWrap, the application first wraps the text to fit the container and then shrinks the text to fit the container.		
	Selections: True or False		
Visible	Specifies whether the selected item appears in the report. By default, the Cover Page and Appendix sections are set to False, and therefore, do not appear in the report.		
	Selections: True or False		

Table 142. Behavior properties in the Properties pane (Sheet 3 of 3)

Data Properties

Table 143 describes the Data properties.

Table 143. Data properties in the Properties pane (Sheet 1 of 3)

Data property	Description		
ColumnCount	(For the Detail Section workspace section only) Specifies the number of columns in the report, similar to a newspaper layout.		
	Default: 1		
	When the template contains data columns that match the layout of the main table (data columns displayed from left to right), use the default value of 1 (ColumnCount = 1) for the number of columns.		
	Tip Increasing the number of columns per page works with transposed data columns (transposed from rows to columns).		
CountNullValues	(For the TextBox design item only)		
	False—Does not include null values as zeros in the summary fields.		
	True—Includes null values as zeros in the summary fields.		
ColumnSpacing	(For the Detail Section workspace section only) Specifies the space between columns (newspaper style) in a multi-column report.		
Culture	(For the TextBox design item only) Formats the report data based on the selected culture from a particular country or region.		
	Default: Inherit		

Data property	Description	
DataField	Specifies the data source (main table column, related table column, or data graph) for the content.	
	IMPORTANT For best results, do not manually modify the content of this box; that is, do not select from the associated dropdown list or type text in the box.	
	To change the design item to another design item, delete the current item. Then, add a new design item as described in "To add a main table column to the report template" on page 455 or "To add a data graph that is associated with the main table to the template" on page 456.	
Description	(For the Picture design item only) Not implemented.	
Hyperlink	Sets to a URL address for a specific location. The application automatically converts this URL to a hyperlink in the HTML or PDF exported reports.	
Image	(For the Picture design item only) Opens the Open dialog box where you can find and select the image file. The default file types for the search are image files.	
MaxLength	(For the RichTextBox design item only) Specifies the maximum number of characters to be displayed.	
NewColumn	(For the Detail Section workspace section only) Specifies where a new column is printed. The default number of report columns is 1. The standard report templates contain only one column and the Customize Reports dialog box creates reports with one column. In this context, a column is a formatting option, not a table column from a result table.	

Table 143. Data properties in the Properties pane (Sheet 2 of 3)

Data property	Description		
NewPage	Specifies whether a page break is inserted before, after, or both before and after the section.		
	Default settings: • CoverPage—After • DetailSection—None • Appendix—Before		
	 CoverPage section: None—No page break between the cover page and the next section. Before—No effect. After—Adds a page break between the cover page and the next section. BeforeAfter—Adds a page break between the cover page and the next section. 		
	 DetailSection: None—No page break between the table rows. Before—Adds a page break between each table row. After—Adds a page break between each table row. BeforeAfter—Adds a page break between each table row. 		
	Appendix section:		
	 None—No page break. Before—Adds a page break before the Appendix section. After—No effect. BeforeAfter—Adds a page break before the Appendix section. 		
Tag	Displays information associated with an object on the page.		
	IMPORTANT Do not modify or delete the Tag property.		
Text	For a Label design item, you can type text here. For the column headings, this box displays the column heading text. For the TextBox design item, this box displays the name of the data source (table column in the result file).		

Table 143. Data properties in the Properties pane (Sheet 3 of 3)

Design Properties

Table 144 describes the Design property.

Table 144.Design properties in the Properties pane

Design property	Description
(Name)	Displays the internal name of an object on the report template page, used by the application to uniquely identify each individual object.

Layout Properties

Table 145 describes the Layout properties.

Table 145.Layout properties in the Properties pane (Sheet 1 of 2)

Layout property	Description	
End	(For the CrossSectionLine and CrossSectionBox design items only) Specifies the X and Y coordinates of the end of the line or the bottom right corner of the box, based on the rulers at the top and left side of the report template page.	
Height	(For a workspace section only) Specifies the height of the section, based on the ruler to the left of the template workspace.	
Location	Specifies the X and Y coordinates of the upper left corner of an object, based on the rulers at the top and to the left of the template workspace.	
Padding	Specifies the values in points for the space to the left, top, right, and bottom of the textual content within the container.	
Radius	(For the CrossSectionBox design item only) Specifies the percentage value for the roundness of the corners of the box. The default value of 0 creates corners with no rounding. A value of 100 creates top and bottom sides that look like half circles.	
RoundingRadius	(For the Shape design item only) Specifies the percentage value for the roundness of the corners when you select RoundRect (rectangle with rounded corners) for the Style property.	
Size	Specifies the width and height of an object on the page, in inches.	
SizeMode	(For the Picture design item only) Specifies how the report designer sizes the image to fit in the container.	
	 Selections: Clip—Clips images that are larger than the container. Stretch—Stretches images to fit the container. Zoom—Decreases the image size to fit the container. 	

Layout property	Description		
Start	(For the CrossSectionLine and CrossSectionBox design items only) Specifies the X and Y coordinates of the start of the line or the top left corner of the box, based on the rulers at the top and to the left of the template workspace.		
X1	(For the Line design item only) Specifies the coordinate of the left end of a line, based on the horizontal ruler at the top of the report template page.		
X2	(For the Line design item only) Specifies the coordinate of the right end of a line, based on the horizontal ruler at the top of the report template page.		
Y1	(For the Line design item only) Specifies the coordinate of the left end of a line, based on the vertical ruler to the left of the template workspace.		
Y2	(For the Line design item only) Specifies the coordinate of the right end of a line, based on the vertical ruler to the left of the template workspace.		

Table 145.Layout properties in the Properties pane (Sheet 2 of 2)

Miscellaneous Properties

Table 146 describes the Miscellaneous properties.

Miscellaneous property	Description	
Code128	(For the Barcode design item only) When you select Code_128_x for the Style property in the Appearance area, use this property to define the code.	
Code49	(For the Barcode design item only) When you select Code49 for the Style property in the Appearance area, use this property to define the code.	
DataMatrix	(For the Barcode design item only) When you select the DataMatrix option for the Style property in the Appearance area, use this property to define the code.	
PDF417	(For the Barcode design item only) When you select Pdf417 for the Style property in the Appearance area, use this property to define the code.	

 Table 146. Miscellaneous properties in the Properties pane (Sheet 1 of 2)

Miscellaneous property	Description	
QRCode	(For the Barcode design item only) When you select QRCode for the Style property in the Appearance area, use this property to define the code.	
	A QR code (quick response code) is a matrix (2D) barcode. Because smartphones can convert QR codes to URLs, QR codes can provide quick access to websites.	
RssExpandedStacked	(For the Barcode design item only) When you select RssExpandedStacked for the Style property in the Appearance area, use this property to define the code.	

Table 146. Miscellaneous properties in the Properties pane (Sheet 2 of 2)

Opening the Property Dialog Boxes for the Report Template's Design Items

The reporting feature for the application includes a property dialog box for each design item in the Section Reports pane. Using this dialog box, you can modify the formatting parameters for a selected design item on the report template page. The available properties vary depending on the selected workspace section or design item. Most of these parameters are similar to the properties listed in the Properties pane, although some might have slightly different names.

* To open the property dialog box for a report design item

1. Select the design item of interest in the workspace.

The properties pane for the selected item appears (Figure 134).

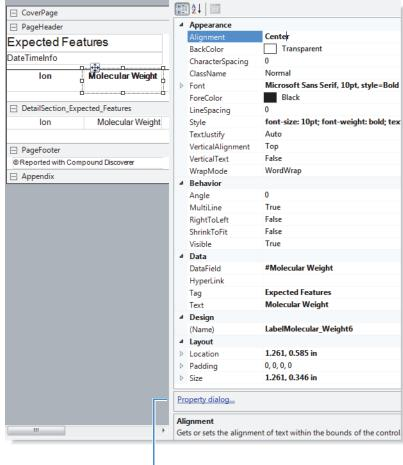


Figure 134. Properties pane for a label design item

Property Dialog link

2. At the bottom right corner of the report template page, click the **Property Dialog** link.

The property dialog box for the selected object opens (Figure 135).

Figure 135. Property dialog box for the Label design item

Label - General	
Label - General General Appearance A Font Format A Alignment	Name: LabelMolecular_Weight6 Tag: Expected Features Visible DataField: #Molecular Weight
	Text: Molecular Weight HyperLink:
	OK Cancel

Note For the RichTextBox design item, in addition to the Property Dialog link, you can click the Load File link to load text from a file into the box (see "Populating a Rich Text Box by Loading a File" on page 487).

The parameters in the property dialog boxes have equivalent parameters in the properties pane.

Populating a Rich Text Box by Loading a File

Use the Load File link at the bottom right corner of the report template page, to populate a RichTextBox design item with data from a file.

To populate a RichTextBox design item

- 1. Click the **Load File** link.
- 2. In the Open dialog box, select the data file and then click **Open**.

You can load text from one of these file types: RTF, TXT, or HTML or HTM.

Selecting the Paper Type, Print Width, Page Orientation, and Watermark for a Report Template

On the report template page, you can add a watermark, change the print width and the orientation of the report template (portrait or landscape), and select the type of paper that you want the printer to use.

Follow these procedures as needed:

- To open the property settings for the report template
- To add a watermark to the report template
- To change the print width of the template
- To change the page orientation of the report template
- To change the paper size used by the printer

Selecting the Paper Type, Print Width, Page Orientation, and Watermark for a Report Template

To open the property settings for the report template *

Click the square in the upper left corner (below the toolbar) of the report template page.

Click this	Target Omeprazole × S Expected Compounds per File ×		
square.			a 🗈 X : 🔉 🖗 100% ▪ 🕸
			. 2
		🖂 CoverPage	
	=	PageHeader	
	-	Expected Compound	ls per File
	<u> </u>	DateTimeInfo	
	=	DetailSection_Expected_Complexity	pounds_per_File
	-	Parent Compound	Parent Compound
	-	Formula	Formula

The watermark properties appear in the Appearance group, and the orientation and paper size properties appear in the Miscellaneous group (Figure 136).

⊿	Appearance							
	ShowParameterUI	True						
	Watermark	(none)						
	WatermarkAlignment	Center						
	WatermarkPrintOnPages							
	WatermarkSizeMode	Clip						
4	Behavior							
	MaxPages	100						
	PrintWidth	8.036						
	ScriptLanguage	C#						
⊿	Data							
	Culture	(default, inherit)						
	DataMember							
\triangleright	DataSource	Thermo.CD.ReportingModule.DataSources.TypedListDataSource						
	UserData	Expected Compounds per File						
⊿	Design							
	TrayHeight	80						
	TrayLargeIcon	False						
⊿	Misc							
	ExpressionErrorMessage							
	IsLandscape	False						
	PaperKind	Α4						
	Version	7.3.7973.0						

Figure 136. Report template properties (in the Properties pane)

Adds a specified image to the report's background. The watermark image can be positioned, sized, aligned and placed on specified pages by using the other watermark properties.

* To add a watermark to the report template

1. Under Appearance, click the browse icon,, to the far right of the Watermark property.

The Open dialog box opens with the file type setting for all image files.

- 2. Find and select the image file of interest and click **Open**.
- 3. In the WatermarkAlignment list, select the appropriate alignment.
- 4. In the WatermarkPrintOnPages box, type the page number of the page where you want the watermark to appear.
- 5. In the WatermarkSizeMode list, select whether you want to clip the image if it is larger than the container, stretch the image to fit the container, or reduce the image size to fit the container.

* To change the print width of the template

Under Behavior, type a numeric value in the PrintWidth box.

* To change the page orientation of the report template

Under Misc, in the IsLandscape list, select True for Landscape or False for Portrait.

* To change the paper size used by the printer

Under Misc, in the PaperKind list, select the appropriate paper size.

To open the list to the paper size of interest, type the first letter of the paper-size name. For example, if you want to change the size from A4 to Letter, type an L in the box. If the paper size does not appear in the box, continue typing the next letter of its name. The application cycles through the paper sizes that begin with this letter or letters.

Previewing and Printing a Report

This topic describes how to work with the Report Preview and Report Print dialog boxes and the report resolution page.

You can print all of the report pages from the report resolution page or the Report Print dialog box. You can print up to five pages from the Report Preview dialog box.

Follow these procedures as needed:

- To open the Report Preview dialog box
- To open the Report Print dialog box
- To open the report resolution page
- To find a text item in a resolved report

- To copy a portion of a report to the Clipboard
- To export the contents of a report to an external document
- To print a report

To open the Report Preview dialog box

In the toolbar on the report template page, click the **Preview Report** icon, **D**.

The application opens the Report Preview dialog box and resolves the current result file with the current template on the report template page. The Page Thumbnails pane displays up to five resolved pages. A red line in the right page margin indicates that the print width does not fit on the selected page size, and the printer will print an empty page with every report page it prints, unless you resize the template (see To change the print width of the template).

Figure 137. Report Preview dialog box with a red line indicating an insufficient right margin

• 弗 오 유, 100 % bnails	• 🕽 🔂 💼 🕅		1/5	J 🖲 🕲	🖑 🕨 🛅			
	Compoun 29-Nov-2015 9	ds per File						Ó
	Molecu	lar Weight	RT [min]		# Adducts	Area		
		345.11458	9.539		1	19124935		
	Urine_0-4hr MW: 345.11	_02, #2404, RT=9. 458, Area: 191249	536 min, FTMS (+) 35					
	(9v0) [stu 250 200 200		346.12186 [M+H]+1					
2	150 [court 100 100 100 100 100 100 100 100 100 100		346.51556	347.12421	348.12726 348.11700	349.12997 349.12033	350.12332	
	10 4	345	346	347	348 m/z	349	350	
3 	(9v0t) [strinoc] viscontra	108.08073 123.0 106.06511 421.08		198	.05827 199.08133 26	7.13641 298.11899 ³¹⁴ 79.10910 312.159419	4.14502 348.12 313.14175	<u> </u>
4	50	100 0.4hz	150		200 m/z	250 3	00	350
	(9,01) (10,00) 300 (10,00) (10,00) 100 (10,00) (10,00) 100 (10,00) (10,00) 100 (10,00)	— 0-4hr	9.536					
	- 0 -	9.4	9.5	9.0 5	3 3 3T [min]	9.7	9.8	
					. found			

To open the Report Print dialog box

In the toolbar on the report template page, click the **Print Report** icon,

The application opens the Report Print dialog box and resolves the current result file with the current template on the report template page. The Page Thumbnails pane displays all of the report pages.

✤ To open the report resolution page

- 1. Open a result file.
- 2. Select an existing report template as follows:
 - a. Choose **Reporting > Create Report** from the menu bar or, click the **Create Report** icon,

The Open Report Design Template dialog box opens to the Report Templates folder.

b. Select the appropriate report template and click **Open**.

The application resolves the current result file with the selected template. The Page Thumbnails pane displays all of the report pages.

To find a text item in a resolved report

- 1. On the report template page, resolve the report by clicking the **Print Report** icon, , or the **Preview Report** icon, .
- 2. In the toolbar of the Report Print dialog box or the Report Preview dialog box, click the **Find** icon, **P**.

The Find dialog box opens.

3. In the Find What box, type the text that you want to find (see Figure 138).

Figure 138. Find dialog box with a text entry

Find	×
Find what: oxidation	Find Next
- w	Find Prev
Match whole word only Match case	Cancel

- 4. Select the appropriate check boxes.
 - To match whole words only, select the Match Whole Word Only check box.
 - To match the case—uppercase or lowercase, select the Match Case check box.
- 5. Click **Find Next** or **Find Prev**.

***** To copy a portion of a report to the Clipboard

Do one of the following:

- a. To copy a specific item to the Clipboard, in the Report Preview toolbar, click the **Selection Mode** icon, **I>**.
- b. Click the item that you want to copy, and then click the **Copy** icon, 💼.

-or-

- a. To copy a rectangular portion of the report to the Clipboard, click the **Snapshot Mode** icon, **1**.
- b. Drag the cursor across the area of interest.

* To export the contents of a report to an external document

- 1. Open the report resolution page or the Report Print dialog box.
- 2. In the toolbar, choose **Export** > *File Type*, where the *File Type* is one of the following:
 - Text—text file
 - PDF—portable document format file
 - RTF—rich text format file
 - Excel—Microsoft spreadsheet
 - HTML—web page that opens in a browser

To print a report

- 1. Do one of the following:
 - a. From the menu bar, choose **Reporting > Create Report**.

The Open Report Design Template dialog box opens.

b. Select an appropriate template and click **Open**.

The report resolution page opens.

-or-

- From the report template page, open the Report Print dialog box by clicking the **Print Report** icon,
- 2. In the toolbar, click the **Print** icon, 🔿 .

The Print dialog box opens.

- 3. Select the appropriate printer and the page range that you want to print.
- 4. Click **OK** to print the report.

Related Topics

- Report Preview Toolbar
- Page Thumbnails Pane

Report Preview Toolbar

The Report Preview and Report Print dialog boxes and the report resolution page share a common toolbar. Table 147 describes the icons in the toolbar, from left to right.

Table 147.	Report	preview	icons	(Sheet 1	of 2)
		p		10001 .	<u> </u>

lcon	Description
R	Toggle Sidebar—Opens and closes the Page Thumbnails pane (see "Page Thumbnails Pane" on page 495).
<u>ا</u>	Print—Opens the Print dialog box where you select the appropriate print options and send the report to the selected printer.
	Copy—Copies the selected item to the Clipboard. Clicking the Selection Mode icon activates the Copy icon.
#1	Find—Opens the Find dialog box where you can search for a particular word or phrase in the report.
2	Zoom Out—Reduces the magnification of the report view. The current zoom box displays the magnification.
8	Zoom In—Increases the magnification. The current zoom box displays the magnification.
100 % -	Current zoom—Use to change the on-screen magnification of the report by selecting or typing a percentage from 10 to 800 in this box, and then pressing ENTER.
 €→	Fit Width—Sizes the width of the report to the screen width. The current zoom box displays the magnification.
\$	Fit Page—Sizes the current report page to the screen width and height, while maintaining the aspect ratio. The current zoom box displays the magnification.
	Single Page View—Fits the current report page to the full-screen view and removes the scroll bar. To view the report pages, you must use the First Page, Previous Page, Next Page, and Last Page icons or the Page Thumbnails pane.
	Continuous View—Changes the magnification to 100% and makes the scroll bar available so that you can scroll through the document.
88	Multipage view—Changes the display to the selected multi-page view.
	First Page—Displays the first page of the report.

lcon	Description
	Previous Page—Displays the previous page of the report.
1/5	Current Page—Indicates the current page and the estimated number of pages in the report.
۲	Next Page—Displays the next page of the report.
	Last Page—Displays the last page of the report.
8	Backward—Displays the previously selected page in the report.
	Selecting pages as you browse activates the Backward icon.
۲	Forward—Displays the next selected page in the report.
	Clicking the Backward icon activates the Forward icon.
4	Pan Mode—Use the hand cursor to drag the page on the screen.
۹I	Selection Mode—Use to copy a report item to the Clipboard.
10	Snapshot Mode—Use to copy a rectangular area to the Clipboard.
Export Text PDF RTF Excel HTML	Export—Use to export the contents of the report to an external document of one of these types: Text, PDF, RTF, Excel, or HTML.

Table 147. Report preview icons (Sheet 2 of 2)

Page Thumbnails Pane

The Page Thumbnails pane appears to the left of the page preview in the Report Preview and Report Print dialog boxes and the report resolution page. Each thumbnail represents a page in the report. The Report Preview dialog box resolves up to five pages of data. The Report Print dialog box resolves all of the data. In the Report Print dialog box, click a page thumbnail to jump to that page.

Table 148 describes the icons in the Page Thumbnails pane, from top left to bottom right.

lcon	Description
+	Enlarge—Enlarges the size of the thumbnails.
-	Reduce—Reduces the size of the thumbnails.
ð	Thumbnails Pane—Displays the Page Thumbnails pane.
æ	Search Results—Displays the Search Results pane where you can search for a particular word or phrase in the report. The found instances appear in the list of results.

Table 148. Page Thumbnails pane icons

11 Creating and Printing Reports

Previewing and Printing a Report

12 -

Using the Lists and Libraries Manager

To modify the lists and libraries that are used to process data in the Compound Discoverer application, see the following topics.

Contents

- Library Manager Page
- Modifying the Expected Compounds List
- Modifying the Adducts List
- Modifying the Ion Definitions List
- Modifying the Transformations List
- Modifying the Mass Lists Library
- Modifying the Spectral Libraries List
- Modifying the Metabolika Pathways List
- Modifying the Compound Classes List
- Loading a Structure from a Structure File
- Finding a Structure in the ChemSpider Database
- Using the Structure Drawing Tools or Commands

Library Manager Page

Clicking the Lists & Libraries Manager icon, \mathbf{m} , in the application toolbar opens the Lists & Libraries Manager as a tabbed page. The buttons in the left pane open the individual views, and the buttons across the top of the page perform various tasks.

Not all the task buttons are available for every view. The Edit button is unavailable for the Spectral Libraries view, and the Replace button is visible only for the Mass Lists, Spectral Libraries, and Metabolika Pathways views.

Each view in the Lists & Libraries Manager page includes a table. For information about freezing panes, hiding and showing columns, freezing rows, sorting, and filtering the application tables, see "Working with the Application Tables" on page 579.

Modifying the Expected Compounds List

The initial expected compounds list contains omeprazole (Figure 139), the targeted compound used in the *Compound Discoverer Metabolism Tutorial*.

Note To use a targeted processing workflow that includes the Generate Expected Compounds node, the Create FISh Trace node, or both of these nodes, you must first add the compounds of interest to the Expected Compounds list.

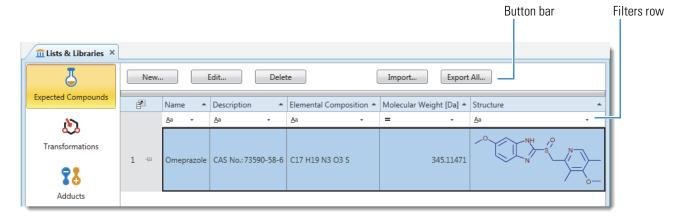


Figure 139. Expected Compounds view

To modify the expected compounds list, see these topics:

- Importing, Exporting, and Deleting Compounds
- Adding and Editing Compounds with the Compound Editor

Importing, Exporting, and Deleting Compounds

Follow these procedures as needed:

- To open the Expected Compounds view
- To delete a compound
- To import a list of compounds from an XML file
- To export the contents of the Expected Compounds view to an XML file

* To open the Expected Compounds view

From the application menu bar, choose **Lists & Libraries > Expected Compounds**.

The Expected Compounds view contains your list of expected compounds for targeted analyses.

To delete a compound

- 1. Select the compound and click **Delete**.
- 2. At the prompt, click **Yes**.

* To import a list of compounds from an XML file

- 1. In the Expected Compounds view, click Import.
- 2. Select the XML file of interest, click **Open**, and then click **OK** at the prompt.

The application only imports new entries; it does not import entries that are already in the list. After the application imports the new compound entries, it provides a tally of the number of imported compounds versus the number of skipped compounds.

* To export the contents of the Expected Compounds view to an XML file

- 1. In the Expected Compounds view, click Export All.
- 2. Select a folder, name the file, and click **Save**.

The application exports the list to an XML file.

3. At the prompt, click **OK**.

Table 149 describes the parameters in the Expected Compounds view.

Table 149. Expected Compounds view parameters (Sheet 1 of 2)

Feature	Description
Buttons	
New	Opens the Compound Editor for adding new compounds to the list.
Edit	Opens the Compound Editor for editing the selected compound.
	Selecting a compound activates this button.
Delete	Deletes the selected compound.
	Selecting a compound activates this button.
Import	Opens the Open dialog box for selecting an XML file.
Export All	Opens the Save As dialog box for exporting the entire list as an XML file.

Feature	Description
Table columns	
Name	Displays the user-specified compound name.
Description	Displays the user-specified description.
Elemental Composition	Displays the elemental composition that the application determines from the compound's structure.
Molecular Weight	Displays the molecular weight that the application calculates from the compound's elemental composition.
Structure	Displays the structure created by using the drawing tools or by importing a structure file.

Table 149. Expected Compounds view parameters (Sheet 2 of 2)

Adding and Editing Compounds with the Compound Editor

Use the Compound Editor dialog box to create new compound entries or edit existing entries in the Expected Compounds view.

Note Similar Compound Editor dialog boxes open from the Mass List and Metabolika Pathways editors. For information about using the structure drawing tools, see "Using the Structure Drawing Tools or Commands" on page 542.

Follow these procedures as needed:

- To open the Compound Editor dialog box
- To add a compound to the Expected Compounds list
- To open a structure file
- To edit an existing entry

To open the Compound Editor dialog box

- 1. Open the Expected Compounds view (see "To open the Expected Compounds view" on page 499).
- 2. Do one of the following:
 - To add a new compound, click New.
 - To edit an existing entry, select the entry and click Edit.

* To add a compound to the Expected Compounds list

1. In the Expected Compounds view, click New.

The Compound Editor dialog box opens . As indicated by the red borders you must name the compound and enter its elemental composition to save it.

Figure 140. Compound Editor dialog box for a new compound

Ocmpound Editor		×
N N N N N / / /	〃∥ r◎○○◎ Დ +-],	
		Click to locate and open a structure file
		Drawing area
Name:		
Description:		
Elemental composition:		7
Molecular weight:		
ChemSpider	Save	
Opens	the ChemSpider application	

2. To enter the elemental composition, draw the structure in the drawing area, open a structure file, or click **ChemSpider** to find the structure.

The application automatically populates the Elemental Composition and Molecular Weight boxes with read-only values. To change the elemental composition and the molecular weight, you must modify the chemical structure in the drawing area.

- 3. (Optional) In the Description box, type a description of the compound.
- 4. Click Save.

The dialog box closes and the compound appears in the list.

✤ To open a structure file

See "Loading a Structure from a Structure File" on page 541.

To edit an existing entry

1. In the Expected Compounds view, either double-click the entry or select it and click Edit.

The Compound Editor dialog box opens with information for the selected entry.

- 2. Edit the information as applicable.
 - To modify the chemical structure, edit the structure in the drawing area (see "Using the Structure Drawing Tools or Commands" on page 542) or open a different structure file.

The application automatically populates the Elemental Composition and Molecular Weight boxes.

- To change the compound name, type alphanumeric text in the Name box.
- 3. To save the changes, click Save, and then click Yes at the prompt.

Modifying the Adducts List

Use the Adducts view to define the adducts in your sample solutions (Figure 141).

Figure 141. Adducts view

îîîLists & Libraries ×							
5	New	Edit		Delete	Impor	t	Export All
Expected Compounds	Ē	Name	▲ A	dduct Mass 🔺	Elemental Composition 🔺	Charge 🔺	
<u>ک</u>		<u>A</u> a •	- =	-	<u>A</u> a •	= -	
~ ~	3 	Ca		39.96149	Ca	2	
Transformations	1 👳	ACN		41.02655	C2 H3 N	0	
	2 🗇	Br		78.91889	Br	-1	
₩	4 ⊹⊐	CI		34.96940	CI	-1	
Adducts	5 👳	DMSO		78.01394	C2 H6 O S	0	
	6 🗇	e		0.00055		-1	
<u>S</u>	7 🗇	FA		46.00548	C H2 O2	0	
Ion Definitions	8 🗇	Fe		55.93330	Fe	3	
1011 Definitions	9 🗇	Н		1.00728	Н	1	
T.	10 🗇	H2O		18.01056	H2 O	0	
Da	11 🕀	HAc		60.02113	C2 H4 O2	0	
Mass Lists	12 🕀	К		38.96316	К	1	
	13 🐤	MeOH			C H4 O	0	
<u>hil.</u> .101	14 👳	Na			Na	1	
Spectral Libraries	15 🗢	NH3		17.02655		0	
-person containes	16 中	NH4		18.03383		1	
●⊥ •	17 🕂	TFA		113.99286	C2 H F3 O2	0	

To modify the adducts list, see these topics:

- Working with the Adducts List
- Adding and Editing Adducts with the Adduct Editor

Working with the Adducts List

Follow these procedures as needed:

- To open the Adducts view
- To delete an adduct from the list
- To import a list of adducts from an XML file
- To export the entire list of adducts to an XML file

* To open the Adducts view

From the application menu bar, choose **Lists & Libraries > Adducts**.

✤ To delete an adduct from the list

- 1. Select the adduct and click **Delete**.
- 2. At the prompt, click Yes.

To import a list of adducts from an XML file

- 1. Click Import.
- 2. In the Open dialog box, locate the file, and click **Open**.

A message opens with a tally of the number of imported adducts versus the number of skipped adducts. The application only imports new entries; it does not import entries that are already in the list.

3. At the prompt, click **OK**.

* To export the entire list of adducts to an XML file

- 1. Click Export All.
- 2. Select a folder, name the file, and click **Save**.
- 3. At the prompt, click **OK**.

Adding and Editing Adducts with the Adduct Editor

Use the Adduct Editor dialog box to define additional adducts, which are part of the ion definition.

To add adducts to the Adducts library

- 1. Open the Adducts view (see "Modifying the Adducts List" on page 502).
- 2. In the button bar, click New.

The Adduct Editor dialog box opens (Figure 142).

Figure 142. Adduct Editor dialog box

Adduct Editor				-
Name:				
New				
Elemental composi	tion:			
Charge:				
0				
		Save	Ca	incel

- 3. Define the new adduct as follows:
 - In the Name box, select **New** and type a name for the adduct.
 - In the Elemental Composition box, type the elemental composition of the adduct. The editor validates the composition entry.
 - In the Charge box, type the charge that the adduct adds to the ion definition.

Range: -10 to +10

4. Click Save.

The new adduct appears in the library.

✤ To edit an adduct entry

- 1. Select the entry.
- 2. Click Edit.

The Adduct Editor opens with the selected entry.

- 3. Modify the adduct definition as follows:
 - In the Name box, select the current name and type a new name for the adduct.
 - In the Elemental Composition box, select the current elemental composition and type a new elemental composition for the adduct.

The editor validates the composition.

• In the Charge box, type the charge that the adduct adds to the ion definition.

Range: -10 to +10

4. Click Save.

Modifying the Ion Definitions List

To modify the Ion Definitions list, see these topics:

- Working with the Ion Definitions List
- Adding or Editing Ion Definitions with the Ion Definition Editor

Working with the Ion Definitions List

The application uses the entries in the Ion Definitions list in the following workflow nodes: Detect Compounds and Generate Expected Compounds.

The default Ion Definitions list contains the most common adducts and dimers formed when using the electrospray-mass spectrometry (ESI-MS) technique in either the positive or negative polarity mode.

Table 150 lists the common adducts and dimers for the positive polarity mode. Table 151 lists the common adducts and dimers for the negative polarity mode.

lon definition	Adducts total mass	Charge	lon definition	Adducts total mass	Charge
M+H–H2O	-17.00329	1	M+H+Na	23.9965	2
M+H–NH3	-16.01927	1	M+H+MeOH	33.03349	1
M+H	1.00728	1	M+K	38.96316	1
2M+H	1.00728	1	2M+K	38.96316	1
M+2H	2.01455	2	M+H+K	39.97044	2
M+3H	3.02183	3	M+H+ACN	42.03383	1
M+NH4	18.03383	1	2M+H+ACN	42.03383	1
2M+NH4	18.03383	1	M+2H+ACN	43.0411	2
M+H+NH4	19.0411	2	M+Na+ACN	64.01577	1
M+Na	22.98922	1	2M+Na+ACN	64.01577	1
2M+Na	22.98922	1	M+H+DMSO	79.02121	1

Table 150. Common adducts and dimers in the positive ion mode

Table 151. Common adducts and dimers in the negative ion mode (Sheet 1 of 2)

lon definition	Adducts total mass	Charge	lon definition	Adducts total mass	Charge
M–H	-1.00728	-1	M–H+FA	44.9982	-1
2М-Н	-1.00728	-1	2M-H+FA	44.9982	-1
M–2H	-2.01455	-2	2M–H+HAc	59.01385	-1

lon definition	Adducts total mass	Charge	lon definition	Adducts total mass	Charge
M-H-H2O	-19.01784	-1	M–H+HAc	59.01385	-1
M+Cl	34.9694	-1	M–H+TFA	112.98559	-1
M–2H+K	36.94861	-1			

 Table 151. Common adducts and dimers in the negative ion mode (Sheet 2 of 2)

Figure 143 shows the default ion definitions.

Figure 143. Ion Definitions view

🟦 Lists & Libraries 🛛 ×						
<u> </u>	New.	Edit Delete	2	Impor	t E	xport All
xpected Compounds	Ē	Ion Definition	Adducts Total Mass 🔺	Charge 🔺	Weight 🔺	
15		<u>A</u> a •		= -	= -	
1	1 👳	2M+H	1.00728	1	50	
Transformations	2 👳	2M+H+ACN	42.03383	1	50	
	3 👳	2M+K	38.96316	1	50	
?	4 ⊹⊐	2M+Na	22.98922	1	50	
Adducts	5 -12	2M+Na+ACN	64.01577	1	50	
Adducts	6 +⊐	2M+NH4	18.03383	1	50	
	7 🖙	2М-Н	-1.00728	-1	50	
N	8 👳	2M-H+FA	44.99820	-1	50	
Ion Definitions	9 +⊨	2M-H+HAc	59.01385	-1	50	
_	10 🗁	M+2H	2.01455	2	50	
	11 👳	M+2H+ACN	43.04110	2	50	
Mass Lists	12 🗇	M+3H	3.02183	3	50	
IVIDSS LISTS	13 🗇	M+CI	34.96940	-1	50	
hill.	14 🗇	M+H	1.00728	1	99	
<u>.dut</u>	15 👳	M+H+ACN	42.03383	1	50	
Spectral Libraries	16 👳	M+H+DMSO	79.02121	1	50	
	17 🕀	M+H+K	39.97044	2	50	
1 1	18 👳	M+H+MeOH	33.03349	1	50	
Metabolika Pathways	19 🗇	M+H+Na	23.99650	2	50	
inclubolika i atiways	20 🗇	M+H+NH4	19.04110	2	50	
• ∩ •	21 🗝	M+H-H2O	-17.00329	1	50	
•~•	22 🗝	M+H-NH3	-16.01927	1	60	
Compound Classes	23 😑	M+K	38.96316	1	60	
	24 😑	M+Na	22.98922	1	70	
	25 🗇	M+Na+ACN	64.01577	1	50	
	26 🖙	M+NH4	18.03383	1	60	
	27 🕂	M-2H	-2.01455	-2	50	
	28 🗇	M-2H+K	36.94861	-1	50	
	29 🗇	M-H	-1.00728	-1	50	
	30 🗇	M-H+FA	44.99820	-1	50	
	31 🕀	M-H+HAc	59.01385	-1	50	
	32 🖙	M-H+TFA	112.98559	-1	50	
	33 中	M-H-H2O	-19.01784	-1	50	

Follow these procedures as needed:

- To open the Ion Definitions view
- To delete an ion definition
- To import a list of ion definitions from an XML file
- To export the contents of the Expected Compounds view to an XML file

To open the lon Definitions view

In the application menu bar, choose **Lists & Libraries > Ion Definitions**.

✤ To delete an ion definition

- 1. In the Ion Definitions view, select the ion definition and click Delete.
- 2. At the prompt, click **Yes**.

To import a list of ion definitions from an XML file

- 1. In the Ion Definitions view, click Import.
- 2. Locate the file and click **Open**.

A message opens with a tally of the number of imported ion definitions versus the number of skipped ion definitions. The application only imports new entries; it skips entries that are already in the list.

3. At the prompt, click OK.

* To export the entire list of ion definitions to an XML file

- 1. In the Ion Definitions view, click Export All.
- 2. Select a folder, name the file, and click **Save**.

The application automatically adds the file name extension (.xml).

3. At the prompt, click OK.

Table 152 describes the columns in the Ion Definitions view.

Table 152. Ion Definitions table columns (Sheet 1 of 2)

Column	Description
Ion Definitions	Displays the user-specified ion definition.
Adducts Total Mass	Displays the difference between the exact mass of the neutral molecule and the molecular ion adduct or the exact mass of the neutral dimer and the ionized dimer.

Column	Description
Charge	Displays the charge of the ion.
Weight	Specifies the weighting factor for the ion definition when the ion definition is added to the list of possible ions in the Ions list for the Detect Compounds node.

Table 152. Ion Definitions table columns (Sheet 2 of 2)

Adding or Editing Ion Definitions with the Ion Definition Editor

Use the Ion Definition Editor dialog box to create new ion definitions.

- ✤ To add ion definitions to the Ion Definitions list
- 1. Open the Ion Definitions view (see "Working with the Ion Definitions List" on page 505).
- 2. In the button bar, click **New**.

The Ion Definition Editor dialog box opens (Figure 144). The Ion Definition box contains only an M for the uncharged molecule. Because the ion definition must include at least one additional component, a charge, or both, the box has a red border, and the Save button is unavailable.

Figure 144. Ion Definition Editor dialog box

Ion Definition Edit	tor 💽
Ion definition:	
М	
Weight factor:	
50	
Available adducts:	
0 📥 ACN	*
0 ≜ Br	≡
0 ▲ Ca	
0 📥 CI	
0 🚔 DMSO	
0 📥 e	
0 🔺 🖬	•
	Save Cancel

3. In the Weight Factor box, type the weighting factor for the ion definition.

Range: 0 to 99

- 4. To enter an ion definition, do any of the following:
 - Type the ion definition in the Ion Definition box.
 - Under Available Adducts, use the arrows to add or subtract components.

-or-

- Select the current value to the left of the component and type a new integer value.
- 5. To save the changes, click Save.

The application calculates the ion's charge and difference in mass between the uncharged molecule and the new ion definition. The new ion definition appears in the list.

To edit an ion definition

- 1. In the Ion Definitions view, select the ion definition.
- 2. Click Edit.

The Ion Definition Editor opens with the selected definition in the Ion Definition box.

- 3. To modify an ion definition, do any of the following:
 - Type the ion definition in the Ion Definition box.
 - Under Available Adducts, use the arrows to add or subtract components.

-or-

- Select the current value to the left of the component and type a new integer value.
- 4. In the Weight Factor box, type a weighting factor from **0** to **99** for the ion.

The Detect Compounds node uses the weight factor value for the ion definitions. With the exception of the protonated molecule $[M+H]^+$ in the positive polarity mode and the deprotonated molecule $[M-H]^-$ in the negative polarity mode, if you set the weight factor to 0, the Detect Compounds node does not look for the specified adduct in the mass spectrum.

5. To save the changes, click **Save**, and then click **Yes** at the prompt.

Table 153 describes the parameters in the Ion Definition Editor dialog box.

Table 153. Ion Definition Editor parameters

Parameter	Description
Ion Definition	Displays the current ion definition. Valid ion definitions include the neutral molecule, which is represented by M, and components from the component list. A red border indicates an invalid ion definition.
	As you edit the ion definition by using the component list, the application automatically updates the ion definition.
Weight Factor	Specifies the weighting factor for the ion definition.
Available Adducts list	Use this list to create custom ion definitions. For information about using this list to create ion definitions, see "To add ion definitions to the Ion Definitions list" on page 508.

Modifying the Transformations List

Follow these topics to modify the Transformations list:

- Deleting, Importing, or Exporting Transformations
- Adding or Editing Transformations with the Transformation Editor

Deleting, Importing, or Exporting Transformations

The Transformations view contains a table of possible transformations (Figure 145). The Generate Expected Compounds node uses a selection of entries from this table and the information in the expected compounds list to generate a table of expected transformations for a known compound.

🟦 Lists & Libraries 🗡									.
5	New.	Edit Del	ete	Import	Export All				
Expected Compounds	₽.	Name *	Leaving Group 🔺	Arriving Group 🔺	Leaving Modification *	Arriving Modification *	ΔM [Da] 🔺	Phase 🔺	Max Occurrence *
2		<u>A</u> a •	<u>A</u> a •	<u>A</u> a •	<u>A</u> a •	<u>A</u> a •	= -	<u>A</u> a	= -
828	1 👳	Acetylation	н	C2 H3 O		C2 H2 O	42.01056	Phase2	1
Transformations	2 🗝	Arginine Conjugation	но	C6 H13 N4 O2		C6 H12 N4 O	156.10111	Phase2	1
	3 🗇	Cysteine Conjugation 1	Н	C3 H6 N O2 S		C3 H5 N O2 S	119.00410	Phase2	1
28	4 ↔	Cysteine Conjugation 2		C3 H7 N O2 S		C3 H7 N O2 S	121.01975	Phase2	1
Adducts	5 👳	Dehydration	H2 O		H2 O		-18.01056	Phase1	2
	6 🕆	Desaturation	H2		H2		-2.01565	Phase1	3
	7 👳	Glucoside Conjugation	н	C6 H11 O5		C6 H10 O5	162.05282	Phase2	1
.0	8 👳	Glucuronide Conjugation	н	C6 H9 O6		C6 H8 O6	176.03209	Phase2	1
Ion Definitions	9 ⊹⊨	Glutamine Conjugation	но	C5 H9 N2 O3		C5 H8 N2 O2	128.05858	Phase2	1
-	10 👳	Glycine Conjugation	но	C2 H4 N O2		C2 H3 N O	57.02146	Phase2	1
Da	11 👳	GSH Conjugation (on Bromine)	Br	C10 H16 N3 O6 S	Br	C10 H16 N3 O6 S	227.15764	Phase2	1
Mass Lists	12 👳	GSH Conjugation (on Chlorine)	CI	C10 H16 N3 O6 S	СІ	C10 H16 N3 O6 S	271.10713	Phase2	1
	13 中	GSH Conjugation (on Fluorine)	F	C10 H16 N3 O6 S	F	C10 H16 N3 O6 S	287.07758	Phase2	1
<u>hili</u>	14 👳	GSH Conjugation 1		C10 H15 N3 O6 S		C10 H15 N3 O6 S	305.06816	Phase2	1
<u>alut</u>	15 🗇	GSH Conjugation 2		C10 H17 N3 O6 S		C10 H17 N3 O6 S	307.08381	Phase2	1
Spectral Libraries	16 中	Hydration		H2 O		H2 O	18.01056	Phase1	1
	17 🗇	Methylation	Н	C H3		C H2	14.01565	Phase2	1
1	18 🗇	Nitro Reduction	O2	H2	02	H2	-29.97418	Phase1	2
Metabolika Pathways	19 😑	Ornitine Conjugation	но	C5 H11 N2 O2		C5 H10 N2 O	114.07931	Phase2	1
inclubolika i dilinayo	20 😑	Oxidation		0		0	15.99491	Phase1	3
0	21 😑	Oxidative Deamination to Alcohol	H2 N	но	HN	0	0.98402	Phase1	1
••••	22 😑	Oxidative Deamination to Ketone	H3 N	0	H3 N	0	-1.03163	Phase1	1
Compound Classes	23 😑	Oxidative Debromination	Br	но	Br	но	-61.91560	Phase1	3
	24 😑	Oxidative Dechlorination	CI	но	CI	но	-17.96611	Phase1	3
	25 😑	Oxidative Defluorination	F	но	F	но	-1.99566	Phase1	3
	26 😑	Palmitoyl Conjugation	н	C16 H31 O		C16 H30 O	238.22967	Phase2	1
	27 😑	Reduction		H2		H2	2.01565	Phase1	1
	28 😑	Reductive Debromination	Br	н	Br	Н	-77.91051	Phase1	3
	29 😑	Reductive Dechlorination	CI	н	CI	н	-33.96103	Phase1	3
	30 🗇	Reductive Defluorination	F	н	F	н	-17.99058	Phase1	3
	31 👳	Stearyl Conjugation	н	C18 H35 O		C18 H34 O	266.26097	Phase2	1
	32 🗇	Sulfation	н	H O3 S		O3 S	79.95681	Phase2	1
	33 😑	Taurine Conjugation	НО	C2 H6 N O3 S		C2 H5 N O2 S	107.00410	Phase2	1
	34 🕀	Thiourea to Urea	S	0	S	0	-15.97716	Phase1	1

Figure 145. Transformations view with the default transformations list

Follow these procedures:

- To open the Transformations view
- To delete an entry from the transformations list
- To import a list of transformations from an XML file
- To export the entire transformations list to an XML file

✤ To open the Transformations view

From the application menu bar, choose Lists & Libraries > Transformations.

***** To delete an entry from the transformations list

- 1. In the Transformations view, select the entry and click **Delete**.
- 2. At the prompt, click Yes.

To import a list of transformations from an XML file

- 1. In the Transformations view, click Import.
- 2. Locate the file and click **Open**.

A message opens with a tally of the number of imported transformations versus the number of skipped transformations. The application only imports new entries; it does not import entries that are already in the library.

3. At the prompt, click OK.

* To export the entire transformations list to an XML file

- 1. In the Transformations view, click Export All.
- 2. Select a folder, name the file, and click **Save**.
- 3. At the prompt, click **OK**.

Table 154 describes the parameters in the Transformations view.

Table 154. Transformations view features (Sheet 1 of 2)

Feature	Description
Buttons	
New	Opens the Transformation Editor dialog box for creating new transformations.
Edit	Selecting an entry activates this button. Opens the Transformation Editor dialog box for editing the selected transformation.

Feature	Description			
Delete	Selecting an entry activates this button.			
	Deletes the selected transformation from the library.			
Import	Opens a dialog box for selecting an XML file to import.			
Export All	Opens the Save As dialog box for saving the list as an XML file.			
Table columns				
Name	Displays the user-specified name for the entry.			
Leaving Group	Displays the user-specified elemental composition of the leaving group for the transformation, if specified.			
Arriving Group	Displays the user-specified elemental composition of the arriving group for the transformation, if specified.			
Leaving Modification	Displays the elemental composition difference between the arriving group and the leaving group, if the transformed compound contains fewer atoms than the original compound.			
Arriving Modification	Displays the elemental composition difference between the arriving group and the leaving group, if the transformed compound contains more atoms than the original compound.			
ΔM [Da]	Displays the difference in mass between the original compound and the transformed compound in daltons.			
Phase	Displays the user-specified category for the transformation.			
Max Occurrence	Displays the user-specified value for the maximum number of times that this transformation can occur in a sequence of combinatorial transformations.			

Table 154. Transformations view features (Sheet 2 of 2)

Adding or Editing Transformations with the Transformation Editor

Use the Transformation Editor to add entries to or to edit entries in the transformations list.

* To add a transformation to the transformations list

1. In the Transformations view, click New.

The Transformation Editor dialog box opens (Figure 146). The empty Name, Arriving Group, and Leaving Group boxes have a red outline. You must enter information in the Name box and in at least one of the group boxes.

Figure 146.	Transformation	Editor	dialog box
-------------	----------------	--------	------------

Transformation Editor	×
Name:	
Arriving group:	
Leaving group:	
Phase:	_
Phase I	•
Maximum occurrence:	
1	
Cancel	2

2. In the Name box, type alphanumeric text to identify the transformation.

The red outline disappears.

- 3. Define the transformation as follows:
 - a. In the Arriving Group box, type alphanumeric text for the arriving group of the transformation, if applicable.

Valid alphabetic characters include all of the naturally occurring elements in the periodic table. The valid range of integers is from 1 to 100 000.

After you define the arriving group, the red outline remains until you place the cursor in the Leaving Group box. If you leave the Arriving Group box empty, the red outline remains until you define the leaving group.

b. In the Leaving Group box, type alphanumeric text for the transformation's leaving group.

The red outline disappears when you select a phase from the Phase list or place the cursor in the Maximum Occurrence box.

- c. In the Phase list, select **Phase 1** or **Phase 2** for a biotransformation, or select **Other** for other transformation types.
- d. In the Maximum Occurrence box, type an integer from 1 to 10.

The Transformation Editor validates the entries from top to bottom.

After you make valid entries, the Save button becomes available.

4. Click Save.

The Transformation Editor dialog box closes and the new entry appears in the transformations library.

* To edit an entry in the transformations list

1. In the Transformations view, select the entry and click **Edit**.

The Transformation Editor dialog box opens. The entry boxes are populated with the information for the transformation that you selected in the transformations list (Figure 147).

Figure 147. Transformation Editor with information for an acetylation chemical reaction

Transformation Editor	×
Name:	
Acetylation	
Arriving group:	
C2 H3 O	
Leaving group:	
Н	
Phase:	
Phase II	-
Maximum occurrence:	
1	
Save	Cancel

2. Make the appropriate changes.

If the changes are invalid, the Save button becomes unavailable, the application outlines the invalid entries in red, and the invalid entries temporarily appear in the transformations library.

3. Click Save and, at the prompt, click Yes.

Table 155 describes the parameters in the Transformation Editor dialog box.

Table 155. Transformation Editor parameters

Parameter	Description
Name	Type alphanumeric text in this box.
Arriving Group	Type the elemental composition of the arriving group in this box.
Leaving Group	Type the elemental composition of the leaving group in this box.
Phase	Select a phase from this list.
	Selections: Phase I, Phase II, and Other
Maximum Occurrence	Type the maximum number of times that this reaction can occur in a set of combinatorial reactions.
	Default: 1 Range: 1 to 10

Modifying the Mass Lists Library

Use the Mass Lists view to create a library of available mass lists to use with the Search Mass Lists node—a processing workflow node that searches selected mass lists for matching compounds. Figure 148 shows the mass lists that come with the application.

To edit the mass lists library or a mass list, follow these topics:

- Adding, Deleting, Importing, Exporting, and Replacing Mass Lists
- Editing a Mass List
- Adding and Editing Mass List Compounds with the Compound Editor

Figure 148. Mass Lists view

Ocmpound Discoverer 3.	0.0.276						
<u>File</u> <u>Reporting</u> <u>Lists</u> & Li	ibraries <u>V</u> iew <u>W</u> indow <u>H</u> elp						
ំ រ៉ារ 🛅 🛅 🕄	🕋 🛱 🏛 🗄 🧃 🌳 : 🞑 🗄 : 🖾 🛗 : 🖳 🗎	V : E 🛛	1 i 🖉	14 : 2, 2, 2, 2	S 6ª 🗎 🕯	Ś.	
🟦 Lists & Libraries 🛛 🗙							~ ×
5	New Edit Delete		Import	Export	Replace		
Expected Compounds	Filename	Description	File Size	Uploaded	Updated	Context	State
2	<u>A</u> a •	<u>A</u> a •	= -	= -	= -	<u>A</u> a •	<u>A</u> a •
	Arita Lab 6549 Flavonoid Structure Database		4,168 KB	5/22/2018 2:54 PM	5/19/2018 12:27 AM		Available
Transformations	EFS HRAM Compound Database		288 KB	5/17/2018 5:27 PM	9/26/2017 4:00 PM		Available
28	Endogenous Metabolites database 4400 compounds		492 KB	5/17/2018 5:27 PM	9/26/2017 3:59 PM		Available
	Extractables and Leachables HRAM Compound Database		1,120 KB	5/17/2018 5:27 PM	9/26/2017 3:59 PM		Available
Adducts							
							
A							
Ion Definitions							
Da Mass Lists							
Da							

The application comes with four mass lists:

- Anita Lab Flavenoid Structure Database.massList (with structures)
- Endogenous Metabolites database 4400 compounds.massList
- EFS HRAM Compound Database.massList
- Extractables and Leachables HRAM Compound Database.massList (with structures)

The application can read the following file types:

- CSV files with a molecular weight column, an elemental composition column, or both of these columns (For best results, verify that the molecular weight is calculated to five decimal places.)
- CSV files created by exporting a ChemSpider Results table or a Compounds table as plain text.

Note The Export > As Plain Text shortcut menu command saves the data in the result table to a CSV file (ChemSpider Results.csv or Compounds.csv).

- CSV files created by exporting data from the Thermo Scientific ToxID application
- CSV files (MaConDa_*version#*_.csv) downloaded from the Mass Spectrometry Contaminant Database
- Mass lists files (.massList) created by the Compound Discoverer application
- XML files created by exporting all the compounds (Export All command) in the Expected Compounds library.

Figure 149 shows a flowchart for adding compounds to or editing the compounds in a mass list.

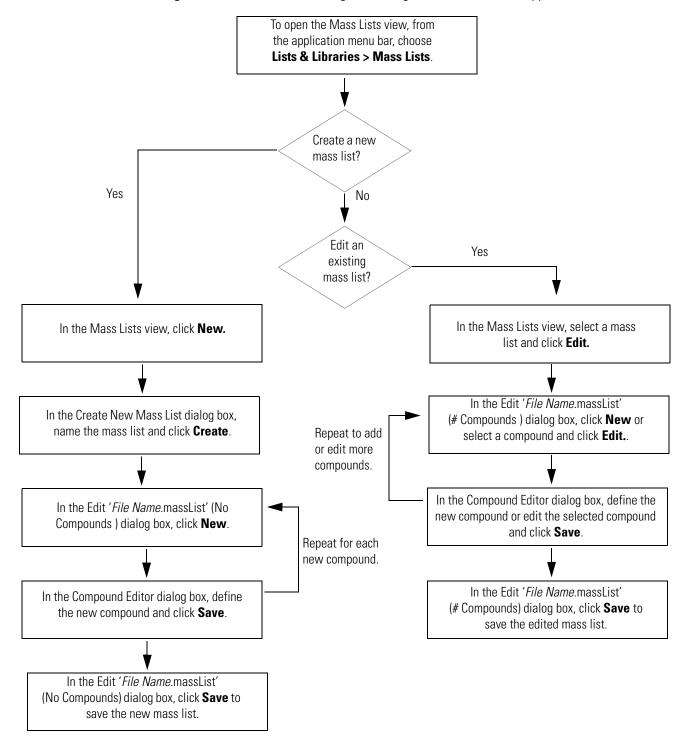


Figure 149. Flowchart for creating and editing mass lists within the application

Adding, Deleting, Importing, Exporting, and Replacing Mass Lists

From the Mass Lists view, you can add, delete, export, and replace mass lists.

Follow these procedures as needed:

- To open the Mass Lists view
- To create a new mass list
- To delete mass list files
- To import a mass list from a CSV file
- To replace a mass list file
- To export a mass list file to a specified file directory

✤ To open the Mass Lists view

From the application menu bar, choose Lists & Libraries > Mass Lists.

✤ To create a new mass list

1. In the Mass Lists view, click New.

The Create New Mass List dialog box opens.

Oreate N	lew Mass List			×
Name:	Required			
Context:				
	<u>(</u>	Create	Cance	:

2. Name the file and click **Create**.

The Edit File Name (No Compounds) dialog box opens.

3. Go to the next topic, "Editing a Mass List" on page 521.

To delete mass list files

- 1. In the Mass Lists view, select the entries to delete and click Delete.
- 2. At the prompt, click Yes.

To import a mass list from a CSV file

- 1. In the Mass Lists view, click Import.
- 2. Locate the CSV file and click **Open**.

The Define CSV File Format dialog box opens.

Note Typically, when the column header in the CSV file matches the name of an essential column, the application recognizes it in the Essential Columns area. A red border around a column name indicates that you must define the column.

The OK button remains unavailable until you define the Name column and at least one column that provides the compound's mass.

- 3. In the Essential Columns area, do the following as needed:
 - To define the column in the CSV file that contains the compound names, select the column name from the Name list.
 - To define the column in the CSV file that provides the compound masses, select the column name from the Formula, Molecular Weight, or Structure lists.

Note Double is the value type for the Molecular Weight column.

- To define the column in the CSV file that provides the chromatographic retention times of the compounds, select the column name from the Retention Time list.
- 4. Click OK.

The Edit 'File Name.massList (# Compounds)' dialog box opens.

5. To edit the mass list, see "Editing a Mass List" on page 521.

To replace a mass list file

- 1. In the Mass Lists view, select the entry and click Replace.
- 2. Browse to the appropriate folder, select a file, and click **Open**.
 - When you select a result file to replace the existing file, the application automatically reads the mass list.
 - When you select a CSV file, the Define CSV File Format for *file name*.csv dialog box opens if the application does not recognize the file format. Otherwise, the application automatically reads the mass list.

To export a mass list file to a specified file directory

- 1. In the Mass Lists view, select the file.
- 2. Click Export.
- 3. Select a folder for the file, rename the file if applicable, and click Save.

Table 156 describes the parameters in the Mass Lists view.

Parameter	Description	
Button		
New	Opens the Create Mass List dialog box for creating and naming a new and empty mass list file. After you name the file and click Create, the Edit ' <i>File Name</i> .massList (# Compounds)' dialog box opens for adding compounds to the list.	
Edit	Opens the Edit ' <i>File Name</i> .massList (# Compounds)' dialog box for editing the selected mass list.	
	Selecting a table row activates this command.	
Delete	Deletes the selected file from the library.	
	Selecting a table row activates this command.	
Import	Opens a dialog box for locating and opening a CSV, a massList, or an XML file that contains, at a minimum, a list of masses.	
	When you select a CSV file, the Define CSV File Format 'File Name.csv' dialog box opens.	
Export	Opens the Save As dialog box for renaming and saving the selected mass list file to another folder.	
Replace	Replaces the selected file with the replacement file. Use this command when the replacement file has the same name as the current file.	
	Selecting a table row activates this command.	
Table columns		
Filename	Displays the file name of the imported file.	
Description	User-editable field for adding descriptive information about the mass list.	
File Size	Displays the file size of the imported file.	
Uploaded	Displays the date (month/day/year) and time (hour/minute) when you added the file to the library in the following format:	
	MM/DD/yyyy HH:mm	
Updated	Displays the data and time when the file was updated.	

Table 156. Mass Lists view parameters (Sheet 1 of 2)

Parameter	Description
Context	Displays the source of the mass list—for example, Import from CSV or Import from XML.
State	Specifies whether the mass list is available, corrupted, or missing.
	If you remove a mass list from the ServerFiles folder or edit a mass list in the ServerFiles Folder, and then restart the application, the mass list's state changes to Missing or Corrupted, respectively.

Table 156. Mass Lists view parameters (Sheet 2 of 2)

Editing a Mass List

Use the Edit '*File Name*.massList (# Compounds)' dialog box to modify the compounds list in a mass list. You can add entries to or edit the existing entries in the list.

To edit a mass list, follow these procedures as needed:

- To open a mass list for editing
- To add a compound to or edit a compound in a mass list
- To delete a compound from a mass list
- To import compounds from an XML file into a mass list
- To import compounds from a CSV file into a mass list
- To add a new column to a mass list

✤ To open a mass list for editing

- 1. From the application menu bar, choose Lists & Libraries > Mass Lists.
- 2. Do one of the following:
 - Create a new mass list (see "To create a new mass list" on page 518).
 - Select a mass list file and click Edit.

-or-

• Double-click a mass list file.

The Edit '*File Name* (# Compounds)' dialog box opens with editable entries for the selected mass list.

* To add a compound to or edit a compound in a mass list

See "Adding and Editing Mass List Compounds with the Compound Editor" on page 524.

✤ To delete a compound from a mass list

- 1. Open the mass list for editing.
- 2. In the Edit '*File Name* (# Compounds) 'dialog box, select the compound, and then click **Delete**.
- 3. Click Save.

* To import compounds from an XML file into a mass list

- 1. Open the mass list (see "To open a mass list for editing" on page 521).
- 2. In the Edit 'File Name (# Compounds) 'dialog box, click Import.
- 3. Locate the XML file and click **Open**.

Note The XML file must have the same format as an XML file created by exporting all the compounds in the Expected Compounds library to an XML file.

4. Click Save.

* To import compounds from a CSV file into a mass list

- 1. Open the mass list (see "To open a mass list for editing" on page 521).
- 2. In the Edit 'File Name (# Compounds) 'dialog box, click Import.
- 3. Locate the CSV file of interest and click **Open**.

The Import Compounds for list 'File Name' dialog box opens. When the application maps a CSV column to an essential mass list column, it automatically displays the column name in the appropriate list. In addition, it makes the check box for a recognized essential column unavailable in the Additional Columns area and selects the data type for the column entries.

- 4. Check the following:
 - In the Essential Columns area, check which CSV columns the application was able to map to the essential columns for a mass list.
 - In the Additional Columns area, check which CSV columns the application was unable to map.

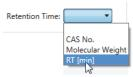
Figure 150 shows the mapped columns from an imported CSV file.

	Mapped columns				
		· · ·			
Import comp	ound for list 'Analgesics'				
Essential Colu	mns				
Name: Nam	e 🔻 Formula: Formula 🔻 Mo	lecular Weight: 📃 🔹 Structure (MOL): Structure Retention Time:		
		, <u> </u>			
Additional Col	umns:				
Use Column	CSV Column Name	Display Name	Value Type		
V	Structure	Structure	String -		
√	Name	Name	String -		
V	Formula	Formula	String -		
V	Molecular Weight	Molecular Weight	String -		
V	RT [min]	RT [min]	String -		
V	CAS No.	CAS No.	String -		
			OK Cancel		
			<u>O</u> K <u>Cancel</u>		

Figure 150. Import Compounds for list 'File Name' dialog box

5. Define any of the remaining columns as appropriate.

For example, to map the RT [min] column in the CSV file as retention time information that the application can search against, in the Essential Columns area, select RT [min] in the Retention Time list.



- 6. (Optional) To avoid importing a non-essential CSV column, clear its check box in the Additional Columns area under Use Columns.
- 7. Click **OK**.
- 8. In the Edit 'File Name (# Compounds)' dialog box, click Save.

To add a new column to a mass list

- 1. Open the mass list (see "To open a mass list for editing" on page 521).
- 2. In the Edit 'File Name (# Compounds)' dialog box, click Add Column.
- 3. In the Add Column dialog box, do the following:
 - a. Name the column.

- b. (Optional) Select the data type:
 - String: Alphanumeric and special characters
 - Double: Decimal numbers
 - Integer: Integers
- c. Click **Add**.
- 4. Click Save.

Table 157 describes the buttons at the top of the Edit '*File Name*.massList' (#Compounds) dialog box.

Table 157. Edit 'File Name.massList' (# Compounds) dialog box parameters

Feature	Description	
Buttons		
New	Opens the Compound Editor dialog box for defining a new compound.	
Edit	Opens the Compound Editor dialog box for editing the selected compound in the mass list.	
	Selecting a compound activates this command.	
Delete	Deletes the selected compound from the mass list.	
	Selecting a compound activates this command.	
Import	Opens a dialog box for locating and opening a CSV, massList, or XML file that contains, at a minimum, a list of masses.	
	When you select a CSV file, the Define CSV File Format 'File Name.csv' dialog box opens.	
Add Column	Opens the Add Column dialog box for naming and defining an additional table column.	

Adding and Editing Mass List Compounds with the Compound Editor

Use the Compound Editor dialog box to add entries to and edit entries in a mass list.

Note The application includes three similar Compound Editor dialog boxes. All three dialog boxes include the same structure drawing tools, but their data entry fields differ. The Compound Editor dialog box that opens for editing mass list compounds includes the following data entry fields: Molecular Weight, Name, Formula, and RT [min].

To add a compound to or edit a compound in a mass list

- 1. From the Mass Lists view, open the mass list for editing (see "Editing a Mass List" on page 521).
- 2. In the Edit 'File Name (# Compounds)' dialog box, do one of the following:
 - To add a compound, click **New**.
 - To edit a compound, double-click it (or select it and click **Edit**).

The Compound Editor dialog box opens (Figure 151). The dialog box is unpopulated for a new compound. The minimum required information is the compound's molecular weight.

Figure 151. Compound Editor dialog box for editing mass list compounds

② Compound Editor	
▶ 🗁 🖬 🗠 🖓 🗸 🖊 🥢	70000 ¢ +,
Molecular Weight:	
0.00000	
Name:	
Formula:	
RT [min]	
	▲ ▼
ChemSpider	Save

3. To change the structure in or add a structure to the drawing area, use the structure tools, load a structure file, or click **ChemSpider** to find a structure file.

For details, see these topics: "Using the Structure Drawing Tools or Commands" on page 542, "Loading a Structure from a Structure File" on page 541, and "Finding a Structure in the ChemSpider Database" on page 541.

Note If you are editing the structure for an existing entry in the mass list, the Molecular Weight field is unavailable, and any formula or structure that you enter must match the displayed molecular weight.

The chemical structure appears in the drawing pane and its molecular weight and formula appear in their respective fields.

4. (Optional) In the Name box, type or edit the compound's name.

5. (Optional) In the RT [min] box, enter a chromatographic retention time.

The application rounds the retention time to three decimal places.

6. Click Save.

The Edit 'File Name (# Compounds)' dialog box appears.

- 7. Review the new or edited compound. If necessary, use the filter row to display only the new entry (see "Filtering the Tables on a Study Page or a List or Library View" on page 588).
- 8. To save the changes to the mass list you are editing, click Save.

Note If you attempt to close the dialog box without saving the changes, an Unsaved Changes prompt appears.

Modifying the Spectral Libraries List

The following mzVault libraries are snapshots of the mzCloud database taken in May 2019:

- mzVault Reference May 2019
- mzVault Autoprocessed May 2019

You can create your own custom mzVault libraries by using the mzVault 2.2 application or by exporting the spectral information from a Compound Discoverer result file (see "Exporting Spectral Data to a New or Existing mzVault Library" on page 286).

Tip The Search mzVault node does not require the mzVault application, so installing the mzVault 2.2 software is optional. However, Thermo Fisher Scientific recommends installing the software to do any of the following:

- Edit existing spectral libraries.
- Create new spectral libraries with curated spectra.
- Convert existing legacy mzVault spectral libraries

You can create new mzVault libraries with the Compound Discoverer 3.1 application, but you are limited to exporting spectra from these result tables: Compounds and Expected Compounds.

Note You cannot edit spectral libraries in the Compound Discoverer application. The Edit button in the command bar is unavailable.

Follow these procedures as needed.

* To open the Spectral Libraries list

From the application menu bar, choose Lists & Libraries > Spectral Libraries.

✤ To add an mzVault library to the list

Click **Import**, locate the library, and click **Open**.

✤ To replace an mzVault library

- 1. In the Spectral Libraries view, select the library.
- 2. Click **Replace**, locate the library, and click **Open**.

To export an mzVault library

- 1. In the Spectral Libraries view, select the library.
- 2. Click **Export**, select a folder, and click **Save**.

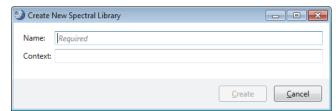
To delete an mzVault library

In the Spectral Libraries view, select the library and click **Delete**.

To create a new empty mzVault library

1. In the Spectral Libraries view, click New.

The Create New Spectral Library dialog box opens.



2. Name the file and click **Create**.

Modifying the Metabolika Pathways List

Initially, the Metabolika Pathways view contains a list of the Metabolika Pathways files (.metabolika) that install with the Compound Discoverer application.

Use the Metabolika Pathways view to create new pathways, edit existing pathways, import and export pathways, and replace pathways.

Follow these topics as needed:

- Deleting, Importing, Exporting, and Replacing Metabolika Pathways
- Editing Metabolika Pathways
- Adding and Editing Pathway Structures with the Compound Editor

Deleting, Importing, Exporting, and Replacing Metabolika Pathways

Follow these procedures as needed.

- To delete pathways
- 1. In the Metabolika Pathways view, select the pathways and click Delete.
- 2. At the prompt, click Yes.

To import pathway files

- 1. In the Metabolika Pathways view, click Import.
- 2. Locate the files (.metabolika) and click **Open**.

To export pathway files to another folder

- 1. In the Metabolika Pathways view, select the pathways, and then click Export.
- 2. In the Browse For Folder dialog box, select a folder and click OK.

The application copies the files to the selected folder.

✤ To replace a pathway

- 1. In the Metabolika Pathways view, select the pathway, and then click Replace.
- 2. Locate the replacement file and click **Open**.

Editing Metabolika Pathways

Use the Edit '*File Name*.metabolika' dialog box to create a new or edit an existing Metabolika pathway.

To create or edit a pathway, follow these topics:

- Creating a New Pathway
- Editing an Existing Pathway
- Modifying Pathway Arrows
- Shortcut Menu for the Metabolika Pathway Editor

Creating a New Pathway

To edit a new (empty) pathway

- 1. From the application menu bar, choose Lists & Libraries > Metabolika Pathways.
- 1. In the Metabolika Pathways view, click New.

The Create New Metabolika Pathway dialog box opens.

2. Name the file and click **Create**.

The Edit Metabolika Pathway 'File Name.metabolika' dialog box opens.

3. In the pathway editor, right-click and choose Add Structure.

Figure 152. Metabolika pathway editor with shortcut menu

🔰 Edit Metabolika Pathway 'New Metabolika Pathway.metabolika' 💼 💷 📧				
Pathway: New Metabolika Pathway.metabolika				
	No data to display	Q 178% A ()		
	Add Arrow			
	Add Structure			
	Add Standard Text	5		
	Merge Text and Arrow	Q ~ [
	Is Reversible			
-	Edit Structure			
	Edit Standard Text			
	Edit Arrow Caption	=		
	Reshape			
	Is Locked			
-	Сору			
	Select All Ctrl+A			
		*		
	III	.]		
		Save		

The Compound Editor opens.

- 4. Use the Compound Editor to add new structures (see "Adding and Editing Pathway Structures with the Compound Editor" on page 533).
- 5. To add an arrow to indicate the direction between reactants and products, right-click a structure and choose **Add Arrow**. To add more arrows, repeat this step.

By default, arrows are straight unidirectional lines with one arrow head and four anchor points (Figure 153). To change the arrow's properties, see "Modifying Pathway Arrows" on page 531.

Figure 153. Selected arrow with four anchor points

- 6. To clean up the drawing, do any of the following:
 - To delete a structure or an arrow, select it and press the Delete key.
 - To delete all the drawing items, right-click and choose **Select All**. Then, press the Delete key.
 - To move a structure or an arrow, select it and hold down the left mouse button to display the move pointer. Then, drag the items to another location.

Editing an Existing Pathway

✤ To edit an existing pathway

1. In the Metabolika Pathways view, select the pathway and click Edit.

The Metabolika pathway editor opens and displays the selected pathway or the beginning section of the pathway in the left pane and the entire pathway at a lower zoom level in the right pane. Use the tools in the upper right of the left pane to change the zoom level, and the selection window in the right pane to display a different section of the pathway in the left pane.

- 2. To modify the pathway, do any of the following:
 - To edit a structure, right-click it and choose Edit Structure.

The Compound Editor opens.

• To add a structure, right-click it anywhere in the left pane and choose Add Structure.

The Compound Editor opens.

- To edit the arrows, see "Modifying Pathway Arrows."
- 3. To undo a change, use the CTRL+Z keys.

You can undo up to six changes.

Modifying Pathway Arrows

To modify the arrows in a Metabolika pathway, follow these procedures as needed:

- To move an arrow
- To change an arrow's angle
- To change an arrow's curvature
- To make an arrow bidirectional to represent a reversible reaction
- To merge arrows
- To add an arrow caption

To move an arrow

1. Select the arrow and hold down the left mouse button to display the move pointer $\binom{+,+}{+}$.



2. Drag the arrow to another location.

To change an arrow's angle

- 1. Select the arrow to display its anchor points, and then point to an end anchor point.
- 2. When this pointer (\mathbf{p}) appears, drag the end of the arrow up or down as needed.

✤ To change an arrow's curvature

- 1. Select the arrow to display its anchor points, and then point to one of its two internal anchor points.
- 2. When this pointer $(\mathbf{\Phi})$ appears, drag the pointer arrow up or down as needed.



✤ To straighten a curved arrow

Right-click it and choose **Reshape**.

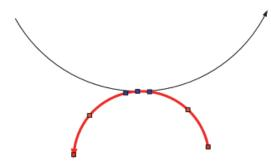
* To make an arrow bidirectional to represent a reversible reaction

Right-click it and choose Is Reversible.

To merge arrows

1. Reshape the arrows into curves.

2. Overlay the curved arrows, and then select one of the curves.



- 3. Point to any of the blue squares.
- 4. When this pointer (**) appears, drag the selected curve until the red square anchor points appear on the second curve. Then, release the mouse button.



5. To straighten one of the merged arrows, right-click it and choose Reshape.

To add an arrow caption

- 1. Right-click the arrow and choose Edit Arrow Caption.
- 2. In the Plain HTML dialog box, type text or text with standard HTML tags and click OK.

Character style	Тад
Bold	
Superscript	
Subscript	

3. To lock the caption to the arrow, right-click the arrow and choose Is Locked.

Shortcut Menu for the Metabolika Pathway Editor

Table 158 describes the shortcut menu for the Metabolika pathway editor.

Table 158.	Shortcut menu	for the	Metabolika	nathway	editor
10010 100.	Unon tout monu		Infotabolika	patrivay	Guitoi

Command	Function
Add Arrow	Adds a straight, unidirectional arrow.
Add Structure	Opens the Compound Editor dialog box for adding a structure.
Add Standard Text	Opens the Plain HTML dialog box for adding formatted text.
Merge Text and Arrow	Merges a selected standard text string to a selected arrow.
Is Reversible	Makes the selected arrow bidirectional.
Edit Structure	Opens the Compound Editor dialog box for editing the selected structure.
	Available for a selected structure.
Edit Standard Text	Use to edit standard text, which is text that is not associated with a structure or an arrow.
Edit Arrow Caption	Opens the Plain HTML dialog box for entering formatted text as an arrow caption.
Reshape	Straightens a curved arrow.
Is Locked	Locks the arrow caption to the arrow.
Copy > Selection	Copies the selected items. Press CTRL+V to paste the copied items elsewhere in the dialog box.
Copy > Structure	Copies the selected structure.
Copy > Pathway	Copies the pathway.
Select All	Selects all the drawing items in the dialog box.

Adding and Editing Pathway Structures with the Compound Editor

Use the Compound Editor to add structures to and edit the structures in a Metabolika pathway.

- ✤ To edit or add a pathway structure
- 1. Open the Compound Editor for editing and adding Metabolika pathway structures (see "Adding and Editing Mass List Compounds with the Compound Editor" on page 524).

Figure 154 shows the Compound Editor dialog box for a new structure.

Ompound Editor	
▶	
	Click to locate and open a structure file.
	Drawing area
Molecular Weight:	
0.00000	
Name:	
Formula:	
ChemSpider	
Opens the ChemSpider application	

Figure 154. Compound Editor dialog box for editing Metabolika pathway structures

- 2. Do one of the following:
 - When adding a new structure, use the drawing tools, load a structure file, or use the ChemSpider application (see "Using the Structure Drawing Tools or Commands" on page 542, "Loading a Structure from a Structure File" on page 541, or "Finding a Structure in the ChemSpider Database" on page 541).

The application automatically populates the Molecular Weight and Formula boxes.

• When editing an existing structure, you are limited to structures that match the current molecular weight and formula.

Modifying the Compound Classes List

When the processing workflow includes the Compound Class Scoring Node, the application can determine the probability that an unknown compound belongs to the user-defined compound classes by comparing the ions detected in the fragmentation scans to user-defined compound classes.

Use the Compound Classes view, the Edit '*Named Compound Class*' dialog box, and the Fragmentation Editor to create a list of compound class libraries. Figure 155 shows the Compound Classes view.

New	Edit Delete		Import	Export		
Name	Description	# Fragments	Uploaded	Updated	File Size	Sta
<u>A</u> a •	<u>A</u> a •	= -	= -		= -	<u>A</u> a
Ecstasy Analogues	MDMA fragments	11	4/25/2017 6:16 PM	4/25/2017 5:28 PM	156 KB	Av
Salicylates	Salicylate fragments	0	4/25/2017 6:02 PM	4/25/2017 6:02 PM	156 KB	Av

Figure 155. Compound Classes view

Follow these procedures as needed:

- To open the Compound Classes library
- To add new compound classes to the Compound Classes library
- To edit a compound class
- To delete a compound class
- To import a compound class library file (.clib)
- To export a compound class

To open the Compound Classes library

From the menu bar, choose Lists & Libraries > Compound Classes.

To add new compound classes to the Compound Classes library

1. In the Compound Classes view, click New.

The Create Compound Class dialog box opens (Figure 156).

Figure 156. Create Compound Class dialog box

Oreate Com	pound Class	
Name:	(Required)	
Description:	(Required)	
		OK Cancel

2. Name the class, provide a brief description, and click **OK**.

The named compound class appears as a new row in the Compound Classes view and the Edit '*Named Compound Class*' dialog box opens (Figure 157). For information about this dialog box, see "Compound Class Editor" on page 538.

Figure 157. Edit 'Named Compound Class' dialog box

🎱 Edit 'Sali	2) Edit 'Salicylates' (No fragments)									
New Edit Delete Import										
r an	m/z		Structure		Formula		Charge	Comment		
	=	•	Aa	*	<u>A</u> a	•	=	<u>A</u> a	•	
									<u>O</u> K	<u>C</u> ancel

- 3. In the Edit 'Named Compound Class' dialog box, click New.
- 4. In the Fragment Editor, define the fragments for the compound class as described in "Using the Fragment Editor" on page 539.
- 5. Click **OK** to close the Edit 'Named Compound Class' dialog box.

To edit a compound class

1. In the Compound Classes view, select the compound class of interest and click Edit.

The Edit 'Named Compound Class' dialog box opens with a list of fragments.

- 2. Do the following as needed:
 - To add a new fragment by using the Fragment Editor, click **New**. Then, use the Fragment Editor to define the fragment (see "Using the Fragment Editor" on page 539).
 - To add new fragments by importing their structures from a CSV or SDF file, click **Import**.
 - To edit a fragment, select it and click Edit.
 - To delete a fragment, select it and click **Delete**.
- 3. Close the dialog box to return to the library.

✤ To delete a compound class

- 1. In the Compound Classes view, select the compound class of interest and click Delete.
- 2. At the prompt, click Yes.

* To import a compound class library file (.clib)

- 1. In the Compounds Class view, click Import.
- 2. Browse to and select a compound class library file (.clib).
- 3. Click **Open**.

The application uploads the file to your local hard drive in the ProgramData > Thermo > Compound Discoverer X.X > Server Files folder. A new entry appears in the library. The entry displays the following information about the compound class: name, description, *#* fragments, date uploaded, date updated, and file size.

To export a compound class

- 1. In the Compounds Class view, select the compound class of interest and click Export.
- 2. In the Save As dialog box, browse to a storage location for the file, change the file name as needed, and click **Save**.

Table 159 describes the features of the Compound Classes view.

Table 159. Compound Classes view parameters (Sheet 1 of 2)

Feature	Description
Commands	
New	Opens the Create Compound Class dialog box for naming and describing the new compound class.
Edit	Opens the Edit ' <i>Named Compound Class</i> ' dialog box for editing the selected compound class file.
Delete	Deletes the selected compound class file.
Import	Opens the Open dialog box for selecting a .clib file.
Export	Opens the Save As dialog box for selecting a file name and location for the exported .clib file.
Table columns	
Name	Displays the name of the .clib file.
Description	Displays a description of the file.
# Fragments	Displays the number of fragments in the file.

Feature	Description
Uploaded	Displays the date (month/day/year) and time (hour/minute) when you added the class to the library in the following format:
	MM/DD/yyyy HH:mm
Updated	Displays the data and time when the file was updated.
File Size	Displays the size of the file.
State	Displays Available when an uncorrupted file with the specified name exists in the following folder:
	<i>Drive X</i> > ProgramData > Thermo > Compound Discoverer <i>X</i> . <i>X</i> > ServerFiles

Table 159. Compound Classes view parameters (Sheet 2 of 2)

Compound Class Editor

Table 160 describes the buttons and columns in the compound class editor. For information about using the compound class editor, see "Modifying the Compound Classes List" on page 535.

Parameters	Description
Buttons	
New	Opens the Fragment Editor for defining a fragment (see "Using the Fragment Editor" on page 539).
Edit	Opens the Fragment Editor with the definition of the selected fragment.
Delete	Deletes the selected fragment.
Import	Opens the Open dialog box for selecting a CSV or an SDF file.
Table columns	
m/z	Displays the m/z value of the fragment.
Structure	Displays the structure of the fragment.
Formula	Displays the formula of the fragment.
Charge	Displays the charge of the fragment.
Comment	Displays the description that you typed in the Create Compound Class dialog box.

Table 160. Edit 'Named Compound Class dialog box parameters

Using the Fragment Editor

Use the Fragment Editor to define fragments for a compound class.

Note For information about using the Fragment Editor's drawing tools, see "Using the Structure Drawing Tools or Commands" on page 542.

To open the Fragment Editor dialog box

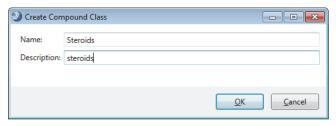
- Open the Compound Classes view from the application menu by choosing Lists & Libraries > Compound Classes.
- 2. In the Compound Classes view, do one of the following:
 - a. Select a library and click **Edit**.

☐ Lists & Libraries ×							~ ×
5	New		Edit	Delete	Import	Export	
Expected Compounds	Name	Description	# Fragments	Uploaded	Updated	File Size	State
0	<u>A</u> a →	<u>A</u> a →	= -	= -		= -	<u>A</u> a →
Compound Classes	Steroids	steroids	0	6/23/2018 3:01 PM	6/23/2018 3:01 PM	156 KB	Available

b. In the Edit 'Named Compound Class' dialog box, select a fragment and click Edit.

-or-

- a. Click New.
- b. In the Create Compound Class dialog box, name the library, type a description, and click **OK**.

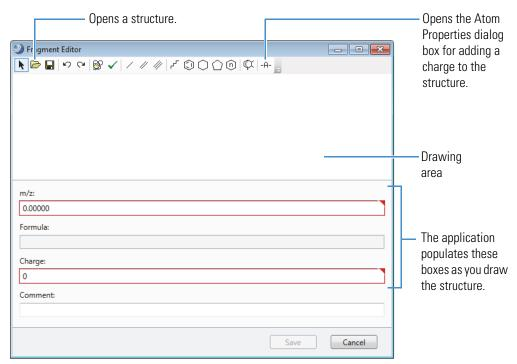


c. In the Edit 'Named Compound Class' (No Fragments) dialog box, click New.

New Edit Delete Import Import Import Import Import <td< th=""><th>② Edit 'Ste</th><th colspan="11"></th></td<>	② Edit 'Ste											
Part Marchard Pormula Charge Comment												
= • <u>A</u> a • <u>A</u> a • = <u>A</u> a •						1			1			
		=	*	<u>A</u> a	*	<u>A</u> a	-	=	<u>A</u> a	-		
	<u> </u>				•		•			•		
									<u>0</u> K	<u>C</u> ance	el	

Figure 158 shows the Fragment Editor dialog box.

Figure 158. Fragment Editor dialog box



* To use the drawing tools to add a structure to the drawing area

See "Using the Structure Drawing Tools or Commands" on page 542.

To open a structure file

See "Loading a Structure from a Structure File" on page 541.

✤ To add a charge to a structure

- 1. In the drawing area of the Fragment Editor, select the atom where you want to add a charge.
- 2. Click the **Atom Properties** icon, ---.
- 3. In the Atom Properties dialog box, select the Charge check box (Figure 159).

Figure 159. Atom Properties dialog box with an available Charge check box

Atom Properties		×
+ C Carbon	Element C H N O F Cl Br I B Si P S R - Substituent	Charge Charge A charge Charge Ch
	Periodic Table	OK Cancel

- 4. Select the positive or negative option.
- 5. Click OK.

Loading a Structure from a Structure File

From the Compound Editor or the Fragment Editor, you can load a structure from a structure file.

- To load a structure from a structure file
- 1. In the editor's toolbar, click the **Load Structure from Disk** button, 🗁.

The Open Structure dialog box opens.

- 2. In the Known Structure Formats list, select the format of the structure file: **MOL Format** (.mol), Compressed Structure (.mcs), or Template (.tml).
- 3. Locate the structure file and click **Open**.

The chemical structure appears in the drawing pane, and the application automatically populates the Elemental Composition and Molecular Weight boxes.

If the structure is not visible or it is only partially visible in the pane, right-click the pane and choose **Select All**. Then, while pressing the SHIFT key, drag the structure into the pane.

Finding a Structure in the ChemSpider Database

You can access the ChemSpider database from the Compound Editor dialog box, the Compound Annotation Editor dialog box, or the mzLogic Analysis view and search for a compound entry with a structure file.

- To load a structure into the drawing area
- 1. Click **ChemSpider**.

The ChemSpider Search dialog box opens.

- 2. In the Input box, enter a name, formula, molecular weight, or CSID.
- 3. Click Search.
- 4. Select a compound from the search results.

Ŋ	ChemSpider Search (Found 1 results)								
In	put: ca	affeine					Search		
CSI	D	Name	Formula	Molecular Weight	# References	Structure			
=	*	<u>A</u> a •	<u>A</u> a -	= -	= -	<u>A</u> a •			
	2424	Caffeine	C8 H10 N4 O2	194.080383	13057				
						Select	Cancel		

5. Click Select.

The chemical structure appears in the drawing pane and the compound's molecular weight, name, and formula appear in their respective fields.

Using the Structure Drawing Tools or Commands

The Compound Editors and the Fragment Editor include a set of drawing tools.

Use the editor toolbar and the shortcut menu for the drawing area to draw, manipulate, and save structures as described in the following topics:

- Compound Editor Toolbar and Fragment Editor Toolbar
- Shortcut Menu Commands for the Drawing Area
- Using the Structure Icons
- Using the Template Tool
- Checking the Validity of a Structure
- Manipulating Structures
- Modifying Atoms and Bonds
- Saving a Structure
- Editing Atom Properties

Note The application has three different Compound Editor dialog boxes and one Fragment Editor dialog box that all have the same structure drawing tools. You can open these dialog boxes from the following Lists & Libraries views: Expected Compounds, Mass Lists, Metabolika Pathways, and Compound Classes.

Compound Editor Toolbar and Fragment Editor Toolbar

The Compound Editor and Fragment Editor dialog boxes have the same set of toolbar icons (Figure 160).

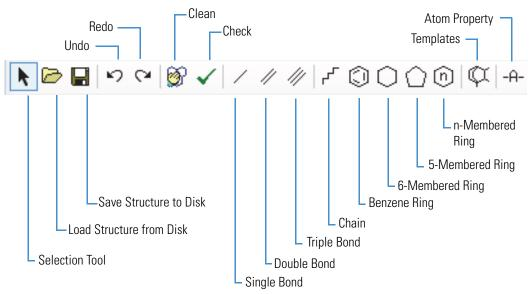


Figure 160. Editor toolbar

Shortcut Menu Commands for the Drawing Area

The Compound Editor dialog boxes and the Fragment Editor dialog box include a drawing area for adding a two-dimensional structure.

Table 161 describes the drawing area's shortcut menu commands.

Menu command	Description	
Selection Tool	Selects a portion of the structure.	
Lasso Selection	Selects an non-rectangular portion of the structure.	
Rectangle Selection	Selects a rectangular portion of the structure.	
Cut	Removes the selected portion of a structure.	
Сору	Copies the selected portion of a structure to the Clipboard.	
Paste	Copies a structure from the Clipboard to the drawing area.	
Delete	Deletes the selected portion of a structure.	
Select All	Selects everything in the drawing area.	
Resize	Resizes the selected portion of a structure (see "To resize a structure" on page 547).	

Menu command	Description
Rotate	Rotates the structure around the selected axis of rotation (see "To rotate a structure" on page 547).
Mirror	Reflects the structure along its vertical or horizontal axis (see "To mirror a structure" on page 547).

Table 161. Shortcut menu commands for the drawing area (Sheet 2 of 2)

Using the Structure Icons

To begin drawing a chemical structure

Tip Point to a drawing icon to display its description.

1. Click any of these structure icons, $/ / / / / / \sim \square \square \square \square \square$.

-or-

Select a template structure as described in "Using the Template Tool."

The cursor changes shape to represent the current drawing mode.

2. Click the drawing area where you want to place the selected structural feature.

Until you click another structure icon, you can continue to add the same structural feature each time you click the drawing area.

3. Edit the atoms and bond properties as described in "Modifying Atoms and Bonds" on page 548.

Using the Template Tool

Use the template tool to draw closely related chemical structures as described in these procedures:

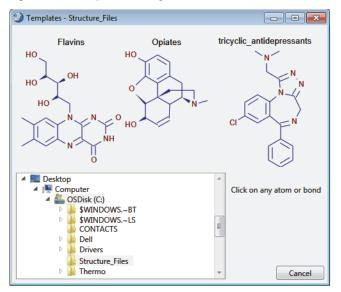
- To open a template structure
- To create a template structure
- To open a template structure
- 1. Click the **Templates** icon, \square .

The Templates dialog box opens.

2. In the Explorer view of the Templates dialog box, browse to and select the folder where you store your structure files.

The title bar of the Templates dialog box changes from Templates to Templates – *Folder name*, and the 2D structures appear above the Explorer view (Figure 161). The application displays all of the structures in the folder. It does not differentiate between MOL files and Template files.

Figure 161. Templates dialog box with a view of the stored structures



3. On the structure that you want to open, click any atom or bond.

The templates cursor, 0, appears in the drawing area of the Compound Editor dialog box.

4. To place the selected structure in the drawing area, click the drawing area.

✤ To create a template structure

- 1. Open a structure file or draw a structure in the drawing area.
- 2. In the Compound Editor toolbar, click the **Save Structure to Disk** button, \square .
- 3. In the Save Structure dialog box, do the following:
 - a. Browse to the folder where you want to store the file.
 - b. Name the file.
 - c. In the Save As Type list, select **Template** (*.tml).
 - d. Click Save.

Checking the Validity of a Structure

The application does not prevent you from creating and saving invalid structures. To check the validity of a structure as you create it or before you save it, use the check structure tool.

Note The check structure tool does not perform quantum mechanical or thermodynamical calculations that address possible structural stability.

✤ To check a structure

1. Click the **Check** icon, \checkmark .

The check structure tool searches for formal errors and unusual structural features. If a structure is formally incorrect or if the check structure tool finds its validity questionable, the Check Structure message lists the errors and warnings (Figure 162).

Figure 162. Check Structure message

Oheck Structure	×
Errors:	
(1) Valence number exceeded on carbon	
Warnings:	
ОК	

2. Click OK.

The application automatically selects the atoms and bonds that it considers incorrect. The application considers structures that are not connected as mixtures and reports them as errors, but it does not select the mixtures.



Manipulating Structures

To manipulate a structure, follow these procedures:

- To resize a structure
- To rotate a structure
- To mirror a structure
- To clean a structure

To resize a structure

- 1. Select the structure or part of the structure you want to resize.
- 2. Right-click and choose Resize.
- 3. Drag one of the small rectangles on the structure's edge and release the mouse button.

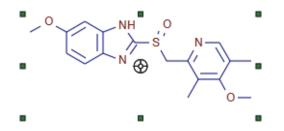
Dragging one of the diagonal rectangles keeps the aspect ratio constant during structure resizing.

✤ To rotate a structure

- 1. Select the structure or part of the structure you want to rotate.
- 2. Right-click and choose Rotate.

A small circle with a cross in the middle, O, appears. The circle indicates the center of rotation.

3. Move the center of rotation by dragging the circle.



4. Rotate the selected structure around the center of rotation by dragging any of the small rectangles on the structure's edge.

To mirror a structure

- 1. Select the structure or part of the structure you want to mirror.
- 2. Right-click and choose **Mirror**.
- 3. Click one of the small rectangles on the structure's edge.
 - The top and bottom rectangles flip the selected structure along a horizontal axis.
 - The left and right rectangles flip the selected structure along a vertical axis.

To clean a structure

- 1. Do one of the following:
 - Select an entire structure.

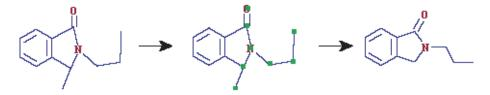
-or-

• Select only the atoms you want to clean.

The selected atoms must be connected.

2. Click the **Clean** icon, \bigotimes .

The cleaning tool helps you create a professional look for your structures.



Note In some complicated cases, the Clean function can lead to structures that you might not find satisfactory. If this occurs, click the **Undo** icon, **S**.

After finishing a structure drawing, always check for errors before proceeding.

Modifying Atoms and Bonds

To modify or add elements to a structure or to modify or add bonds to a structure, follow these procedures:

- Selecting Atoms and Bonds
- Editing Bond Properties

For information about editing the atom properties, see "Editing Atom Properties" on page 550.

Selecting Atoms and Bonds

You can select individual atoms and bonds, a contiguous portion of a structure, an entire structure, or groups of atoms and bonds that are not adjacent to each other.

Follow these procedures:

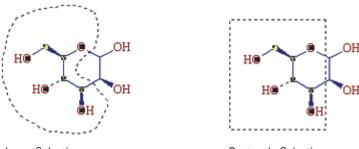
- To select an individual atom or bond
- To choose a selection mode
- To select a group of atoms that are adjacent to each other
- To select all of the atoms and bonds in the structure
- To move a structure

✤ To select an individual atom or bond

Click the Selection Tool icon, and then click the individual atom or bond.

✤ To choose a selection mode

Right-click anywhere in the drawing area of the Compound Editor dialog box and choose **Lasso Selection** or **Rectangle Selection**.



Lasso Selection

Rectangle Selection

* To select a group of atoms that are adjacent to each other

Do one of the following:

• Right-click and choose **Rectangle Selection**. Then drag the cursor to form a rectangle around the atoms.

-or-

• Right-click and choose **Lasso Selection**. Then draw a free-form shape around the atoms.

* To select all of the atoms and bonds in the structure

Do one of the following:

• Right-click the drawing area and choose Select All.

-or-

• Click the **Selection Tool** icon, and then double-click anywhere in the drawing area, except on atoms or bonds.

To move a structure

- 1. Select the atoms or bonds that you want to move.
- 2. Drag the selected structures to a new location.

✤ To move all of the structures in the drawing area

- 1. Right-click the drawing area and choose Select All.
- 2. Click any atom or bond in the drawing area, and then drag the structures to a new location.

Editing Bond Properties

Use the bond icons to change the bond multiplicity.

✤ To change the multiplicity of a bond

Click /, /, or //, and then click the bond that you want to change.

Saving a Structure

After you draw or modify a structure, you can save the structure as a structure file (in MOL format or as a compressed structure) or as a template file.

To save a structure as a structure file

- 1. Click the **Save Structure to Disk** button in the toolbar,
- 2. In the Save Structure dialog box, do the following:
 - a. Browse to the directory where you want to store the file.
 - b. Name the structure file.

You can save structures under their actual names, regardless of length (for example, 1-Amino-2-hydroxyindane.mol).

- c. In the Save As Type list, select a file type.
- d. Click Save.

Editing Atom Properties

Use the Atom Properties dialog box to change the isotope of an atom or the entire element. For more information about modifying compound structures, see "Using the Structure Drawing Tools or Commands" on page 542.

IMPORTANT The application does not support compounds with a charge or radical nor does it support the R-Substituent feature.

* To edit the element or nucleon number of a single atom

- 1. Open the Compound Editor dialog box as described in "To open the Compound Editor dialog box" on page 500.
- 2. Click the **Selection Tool** icon,
- 3. Do one of the following:
 - Double-click the atom that you want to change.

-or-

• Select the atom that you want to change and click the **Atom Properties** icon, -A-.

The Atom Properties dialog box opens with the properties of the selected atom displayed. The Charge and Radical check boxes are not available for compounds.

Tip To add charges to fragment structures for a Compound Class library, see "Using the Fragment Editor" on page 539.

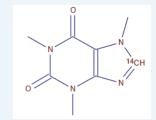
Figure 163. Atom Properties dialog box, showing the properties for carbon

Atom Properties	Element C H N O F Cl Br I B Si P S R - Substituent Periodic Table	Charge Charge C + O - Radical I Isotope Nucleon Number: 12	The Charge and Radical check boxes are not available for compounds.
	I	OK Cancel	

- 4. To change the element, do the following:
 - To change the atom to an element that is in the Element area, click the appropriate Element button.
 - To change the atom to an element not listed in the Element area, click **Periodic Table**. Then, in the Periodic Table dialog box, select an element and click **OK**.
- 5. To specify a less abundant isotope of the element, select the **Isotope** check box, and then select the appropriate value in the Nucleon Number list.

Tip For example, to create a compound that is labeled with one carbon-14 atom, double-click the labeling site—the atom that you want to change. In the Atom Properties dialog box, select the **Isotope** check box, and then select **14** in the Nucleon Number list.

The application displays carbon-14 as [14]C; that is, the elemental composition of carbon-14 labeled caffeine is displayed as C7 [14]C H10 N4 O2.



6. When you finish editing the selected atom, click **OK**.

Changes you make in the Atom Properties dialog box affect only the selected atom.



Using the License Manager

Use the License Manager to activate or deactivate the software license and to install new processing workflow nodes as they become available.

Contents

- Opening the License Manager
- License Manager Command Bar
- Activating the Software License
- Deactivating the Software License for Transfer to Another Computer
- Installing or Updating a Processing Workflow Node

Note After you install the Compound Discoverer application on your computer, you can use the application without activating the license for up to 60 days.

After you activate the software license on one computer, you can deactivate the license and transfer it to another computer.

Opening the License Manager

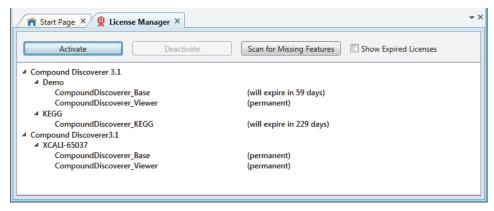
Access the License Manager page from the Help menu.

To open the License Manager page

From the menu bar, choose **Help > License Manager**.

The License Manager opens as a tabbed document in the application window (Figure 164).

Figure 164. License Manager page



License Manager Command Bar

Table 162 describes the License Manager command bar.

Table 162. License Manager command bar

Command or feature	Description
Activate	Opens the License Activation dialog box where you can apply a new activation code and activate the license on the current computer.
Deactivate	Opens the License Deactivation dialog box for deactivating the software license.
Scan for Missing Features	Activates a scan for newly installed processing workflow nodes.
Show Expired Licenses	Selecting this check box displays any expired licenses.

Activating the Software License

To activate your Compound Discoverer 3.1 software license, see these topics as necessary:

- 1. Entering the Product ID and the Activation Code
- 2. Follow one of these topics:
 - Activating the License on an Online Computer
 - Activating the License on an Offline Computer

Entering the Product ID and the Activation Code

To activate your Compound Discoverer 3.1 license, you must know the product ID (XCALI-*XXXXX*) and the activation code. Within one week of ordering the software, you should receive an email from Thermo MS Licensing with the subject line "Your Order Is Ready." This email contains the product ID and activation code information.

✤ To enter the licensing information

1. On the License Manager page, click Activate.

The License Activation dialog box opens to the Activation Code view.

Figure 165. License Activation dialog box opened to the Activation Code view

License Activation	×		
Activation Code			
Locate the email from Thermo MS Licensing with "Your Order is Ready" in the subject line. This email contains the product ID and activation code for your software. If you cannot locate this email, click Help for more information.			
Enter the company nar one of the following:	me, your full name and email address, and the product ID and activation code. Then, do		
 If this computer co 	nnects to the Internet, click Online Activation.		
	bes not connect to the Internet, click Offline Activation and follow the instructions. the response file (activation.xml), click Process Response File on this computer to ation process.		
Company:	Thermo Fisher Scientific		
Full Name:			
User Email:			
Product ID:	kcali		
Activation Code:	<u>·</u> _·_		
Help	Online Activation Offline Activation Process Response File Cancel		

- 2. If you have not already received your activation code, do the following:
 - a. Check your Junk Email folder.
 - b. If the email is not in your Junk Email folder, log in to your account at the following URL. In the left navigation pane, under Software & Services, click **Order History**. Then, in the list of ordered products, click the order number.

https://thermo.flexnetoperations.com

c. If you cannot find your account, send an email message to Licensing at ThermoMSLicensing@thermofisher.com.

Provide the following information in the body of the message:

- Software application: Compound Discoverer
- Sales order number or purchase order number:______
- End user name:_____
- End user email:_____
- 3. In the License Activation dialog box (Figure 165), enter the following:
 - Your company name
 - Your full name
 - Your contact email address
 - The product ID for the Compound Discoverer 3.1 application.

There are five possible product IDs. Four of the product IDs are for software upgrades (see Table 163).

• The activation code. You can type or paste the activation code.

Table 163. Product IDs for the Compound Discoverer 3.1 software (Sheet 1 of 2)

Material Order No.	Product ID	Description
OPTON-30925	XCALI-65037	SW, Compound Discoverer 3.1 (single license)
OPTON-30926	XCALI-65037	SW, Compound Discoverer 3.1 (2–5 licenses)
OPTON-30927	XCALI-65037	SW, Compound Discoverer 3.1 (6–10) licenses
OPTON-30928	XCALI-65040	SW, Compound Discoverer 3.1 upgrade from MetWorks (MW) or Mass Frontier (MF)
OPTON-30929	XCALI-65041	SW, Compound Discoverer 3.1 upgrade from Sieve

Material Order No.	Product ID	Description
OPTON-30930	XCALI-65042	SW, Compound Discoverer (CD) 3.1 upgrade from CD 1.0, CD 2.0, CD 2.1
OPTON-30931	XCALI-65043	SW, Compound Discoverer 3.1 upgrade from Compound Discoverer 3.0

Table 163. Product IDs for the Compound Discoverer 3.1 software (Sheet 2 of 2)

- 4. Depending on whether your processing computer is connected to the Internet, continue with one of these topics:
 - Activating the License on an Online Computer
 - Activating the License on an Offline Computer

Activating the License on an Online Computer

Follow these instructions if your processing computer has an Internet connection.

- * To activate the software license on an online computer
- 1. If you have not already entered the licensing information, enter it in the Activation Code view of the License Activation dialog box.
- 2. Click **Online Activation** to process the activation code.
- 3. At the confirmation prompt, click **OK**.
- 4. At the "new features" prompt, click OK.

This completes the online license activation process.

Activating the License on an Offline Computer

Follow these instructions if your processing computer is not connected to the Internet.

Note Activating the license on an offline computer is a three-step process that requires access to an online computer.

- 1. Create an activation request file (Activation-*Activation Code*.req) on the offline computer.
- 2. Transfer the activation request file to an online computer where you upload it to the licensing portal to obtain a response file (activation.xml).
- 3. Transfer the response file (activation.xml) to the offline computer, and then process it by clicking Process Response File in the License Activation dialog box.

* To activate the software license on an offline computer

- 1. If you have not already entered the licensing information in the Activation Code view of the License Activation dialog box, enter it now.
- 2. In the License Activation dialog box, click **Offline Activation**.

The Save Activation Request File dialog box opens.

3. Save the activation request file (Activation-Activation Code.req).

Save Request File		
😋 🔍 🛡 📕 « Do	ocuments	2
File <u>n</u> ame:	Activation 4444-EEEE-AAAA-5555.req	
Save as <u>t</u> ype:	Request File (*.req)	
Browse Folders	Save Cancel	

The Offline Activation instructions appear.

Figure 166. Offline activation instructions

License Activation	
Offline Activation	
You saved the activation request (.req) file to the following folder: :	
C:\Users\Documents\ACTIVATION REQUEST FILE\Activation 4444-EEEE -AAAA-5555.req	
1. Transfer the activation request file to an online computer, and then go to the following URL:	
https://thermo.flexnetoperations.com/control/thmo/offlineActivation	
2. On the online computer, process the activation request file, and then save the downloaded	
response file (activation.xml).	
3. Transfer the response file to this offline computer, and then click Process Response File.	
Tip: Closing this dialog box copies the URL to the Clipboard.	
	ОК

4. Click OK.

The Activation Code view of the License Activation dialog box reappears.

Tip When you close the Offline Activation dialog box, the application automatically saves the URL to the Clipboard, so that you can save the URL to a file for use on an online computer.

- 5. (Optional) Save the URL to a file.
- 6. Keep the License Activation dialog box open on this offline computer.

IMPORTANT If you accidentally close the License Activation dialog box, start over at step 1.

- 7. To download the response file (activation.xml) from the licensing server, do the following on an online computer:
 - a. Transfer the activation request file ((Activation-*Activation Code*.req) to the online computer.
 - b. Go to the following URL (case sensitive):

https://thermo.flexnetoperations.com/control/thmo/offlineActivation

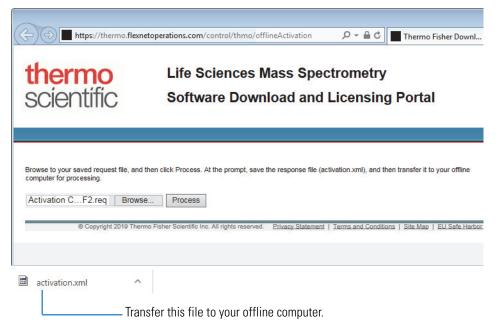
The Life Sciences Mass Spectrometry Software Download and Licensing Portal opens.

c. Click **Choose File** or **Browse**, browse to and select the request file (Activation *Activation Code.*req), click **Open**, and then click **Process**.

Note Whether the page includes a Choose File button or a Browse button depends on the web browser.

The server downloads the activation file.

Figure 167. Software licensing portal with a view of a downloaded response file



d. Save the response file (activation.xml).

- 8. To activate the license on the offline computer, do the following:
 - a. Transfer the response file (activation.xml) to this computer.
 - In the License Activation dialog box that you left open, click Process Response File.
 The Open Response File dialog box opens.
 - c. Browse to and select the response file (activation.xml), and then click Open.
 - d. At the "license response processed" prompt, click OK.
 - e. At the "new features" prompt, click OK.

This completes the offline license activation process. The License Manager indicates that the license for the specified product (XCALI-*XXXXX*) is permanent.

Deactivating the Software License for Transfer to Another Computer

To transfer the software license to another computer, you must first deactivate the license on the current computer.

Follow the appropriate topic:

- Deactivating the Software License on an Online Computer
- Deactivating the Software License on an Offline Computer

Deactivating the Software License on an Online Computer

Follow this procedure if your computer has Internet access.

- * To deactivate the software license on an online computer
- 1. On the License Manager page, select the software product that you want to deactivate, and then click **Deactivate**.

The License Deactivation dialog box opens. The Product ID box is populated with the selected product's ID, and Activation Code box is populated with the activation code for your license.

Figure 168. License Deactivation dialog box populated with the product ID and activation code

Uicense Deactivation	
License Deactiva	ation
To install this license o	n a different computer, you must deactivate the license on this computer.
To deactivate the licen	se, select it, and then do one of the following:
 If this computer co 	nnects to the Internet, click Online Deactivation.
• If this computer do	es not connect to the Internet, click Offline Deactivation. Follow the
instructions to dow	nload a response file on an online computer. Then, process the response file
on this computer.	
Product ID:	XCALI-65037
Activation Code:	4A6C-E7DD-AA4B-575B
Help	Online Deactivation Offline Deactivation Process Response File Cancel

- 2. If the computer is connected to the Internet, click **Online Deactivation**.
- 3. At the confirmation prompt, click Yes.

The license deactivation process is complete, and the product ID disappears from the License Manager page.

Deactivating the Software License on an Offline Computer

If your processing computer is not connected to the Internet, follow these instructions to deactivate the software license.

Note Deactivating the license on an offline computer is a three-step process that requires access to an online computer.

- 1. Create a deactivation request file (Deactivation-*Activation Code*.req) on the offline computer. This step only starts the license deactivation process.
- 2. Transfer the deactivation request file (Deactivation-*Activation Code*.req) to an online computer where you upload it to the licensing portal to obtain a response file (activation.xml).
- 3. Transfer the response file to the offline computer, and then process it to complete the deactivation process.

* To deactivate the software license on an offline computer

- 1. On the offline computer, create the deactivation request file as follows:
 - a. On the License Manager page, select the software license that you want to deactivate, and then click **Deactivate**.

The License Deactivation dialog box opens. The Product ID box is populated with the selected product's ID, and Activation Code box is populated with the activation code for your license.

- b. Click Offline Deactivation.
- c. At the confirm deactivation prompt, click Yes.

The Save Request File dialog box opens. The File Name box displays the following text: Deactivation *Activation Code*.req.

- d. Save the deactivation request file.
- e. At the prompt, which displays the location of the saved request file, click OK.

The license state changes to Deactivation in Progress, and the URL is copied to the Clipboard.

- f. (Optional) Copy the URL to a file.
- 2. From the online computer, retrieve the response file (activation.xml) as follows:
 - a. Transfer the deactivation request file and the file with the URL to this computer.
 - b. Go to the following URL (case sensitive):

https://thermo.flexnetoperations.com/control/thmo/offlineActivation

The Life Sciences Mass Spectrometry Software Download and Licensing Portal opens.

c. Click **Choose File**, browse to and select the **Deactivation** *Activation Code*.req file, click **Open**, and then click **Process**.

The server downloads the response file.

- d. Save the response file (activation.xml).
- 3. Complete the deactivation process on the offline computer as follows:
 - a. Transfer the response file (activation.xml) to this computer.
 - b. On the License Manager page, select the license that you are in the process of deactivating, and click **Deactivate**.

The License Deactivation dialog box opens.

c. Click Process Response File.

The Open Response File dialog box opens.

d. Select the response file (activation.xml) and click Open.

The license disappears from the License Manager page.

Installing or Updating a Processing Workflow Node

The application uses a node-based workflow to process raw data files. Following set guidelines, you can create your own custom workflow nodes. In addition, Thermo Fisher Scientific might occasionally provide custom workflow nodes on its customer website.

* To install a new processing workflow node

- 1. Download the executable files and store them in the appropriate folder on the computer where you are running the application.
- 2. Open the License Manager page.
- 3. Click Scan for Missing Features.
- 4. Close and reopen the application.
- 5. Choose **Help > About**.

The About Compound Discoverer dialog box opens with the Patent and Legal Notices page displayed.

6. Expand the Nodes list and verify that it lists the new node.

13 Using the License Manager Installing or Updating a Processing Workflow Node

Setting Up the Configuration Options

To set up the application's configuration options, follow these topics.

Contents

- Opening the Configuration Page
- Selecting the Maximum Number of Parallel Processing Jobs
- Selecting Where to Store Temporary Data
- Selecting the Global Color Palettes
- Turning Off the Auto-Save Feature for Studies
- Hiding the Workflow Node Numbers
- Specifying the Default mzCloud Mass Tolerance Settings
- Setting Up a BioCyc Account or Subscription
- Specifying the Fragmentation Libraries

Opening the Configuration Page

Use the Configuration page to set up the application's global configuration options.

To open the Configuration page

From the menu bar, choose **Help > Configuration**.

Selecting the Maximum Number of Parallel Processing Jobs

Use the Parallel Options view of the Configuration page to specify the maximum number of analyses (jobs) that the application can process in parallel.

* To change the maximum number of parallel jobs

1. In the left pane of the Configuration page, under Server Settings, select **Parallel Options**.

2. In the Parallel Options view, in the Maximum Number of Processing Workflows in Parallel Execution box, type or select an integer from **1** to **4**.

The default value is equal to half the number of CPU cores in the processing computer.

3. Click Save Current Settings.

Selecting Where to Store Temporary Data

Use the Scratch Directory Options view of the Configuration page to change the folder where the application stores temporary data during data processing.

- ✤ To change the scratch directory
- 1. In the left pane of the Configuration page, under Server Settings, select **Scratch Directory Options**.

The Current Scratch Directory box lists the current location of the scratch folder.

Figure 169. Scratch Directory Options view

Sconfiguration ×			▼ ×
 Server Settings Parallel Options Scratch Directory Options Client Settings (I) Study Management Settings Workflow Editor Settings Color Schema Miscellaneous Settings Submit Single Spectrum to mzCloud Options BioCyc User Login HighChem Fragmentation Options 	Current scratch directory: New scratch directory: Note:	e scratch directory used to store temporary processing data. C:\ProgramData\Thermo\Compound Discoverer 3.0\Scratch	
		🛃 Save Current Setting	gs

- 2. Click the browse button next to New Scratch Directory and locate the new directory.
- 3. Click Save Current Settings.
- 4. Restart the application.
- 5. Reopen the Scratch Directory Options view and make sure that the Current Scratch Directory box lists the new scratch directory.

Turning Off the Auto-Save Feature for Studies

If you want the application to automatically save changes that you make to studies, including the list of result files on the Analysis Results page, do not turn off the auto-save feature.

✤ To turn off the auto-save feature

1. In the left pane of the Configuration page, under Client Settings, select **Study Management Settings** (Figure 170).

Figure 170. Study Management Settings view with the default setting

🛠 Configuration 🗡	• X
 Server Settings Parallel Options Scratch Directory Options Client Settings Study Management Settings Workflow Editor Settings Color Schema Miscellaneous Settings Submit Single Spectrum to mzCloud Options BioCyc User Login HighChem Fragmentation Options 	Study Management Settings Presets affecting the study management Study Related Settings Image: Save study automatically on close and after analysis submit
	Save Current Settings

- 2. Clear the Save Study Automatically On Close and After Analysis Submit check box.
- 3. Click Save Current Settings.
- 4. Restart the application.

Hiding the Workflow Node Numbers

When you create a processing workflow by dragging the workflow nodes into the Workflow Tree pane, the application automatically adds an integer to each workflow node. Use the Workflow Editor Settings view to hide these numbers.

* To hide the workflow node numbers

- 1. In the left pane of the Configuration page, under Client Settings, select **Workflow Editor Settings**.
- 2. Under Workflow Settings, select the Hide Node Numbers check box (Figure 171).



 Server Settings Parallel Options Scratch Directory Options Client Settings (I) Study Management Settings Workflow Editor Settings Color Schema Miscellaneous Settings Submit Single Spectrum to mzCloud Options BioCyc User Login HighChem Fragmentation Options 	Workflow Editor Settings Presets affecting the workflow editor Workflow Settings Image: Image
--	---

3. Click Save Current Settings.

Selecting the Global Color Palettes

Use the Color Schema view of the Configuration page to select one of the standard global color palettes or create a custom color palette. The colormap selection affects the sample group colors in the following views: Chromatograms, Trend Charts, Principal Component Analysis, and Descriptive Statistics. The selection does not affect the color-coding in the result tables.

Figure 172. Color Schema view

	🖉 😽 Configuration 🛛 🕹	
Image: Scratch Directory Options Image: Scratch Directory Options	 Scratch Directory Options Elient Settings Study Management Settings Workflow Editor Settings Color Schema Miscellaneous Settings Submit Single Spectrum to mzCloud Options BioCyc User Login 	Color schema used to colorize data in charts. Selected Palette: Compound Discoverer Keww. Edit Delete Import Export

To set up the global color palette, follow these topics as needed:

- Opening the Color Schema View
- Selecting a Standard Color Palette
- Creating Custom Color Palettes
- Deleting Custom Color Palettes
- Importing Custom Color Palettes

- Exporting Custom Color Palettes
- Editing a Custom Color Palette
- Selecting a Color in the Gradient Color Chart

Opening the Color Schema View

- ✤ To open the Color Schema view
- 1. From the menu bar, choose **Help > Configuration**.
- 2. In the left pane of the Configuration page, under Client Settings, select Color Schema.

Selecting a Standard Color Palette

For visualizing chart data, you can select from four standard color palettes.

To select a standard color palette

- 1. Open the Color Schema view.
- 2. In the Selected Palette list, select from four palettes.
 - Compound Discoverer



• Deuternopia, Protanopia, and Tritanopia

3. Click Save Current Settings.

Creating Custom Color Palettes

- To create a custom color palette
- 1. Open the Color Schema view.
- 2. Click New.

The Color Palette Configuration dialog box opens.

② Color Palette Configuration		
Palette name:		
Palette colors:		
	Red: Green: Blue: Hue: Saturation: Lightness: Hexcode:	255 0 0 0 100% 50% #FF0000
Advanced options	Preview:	
Add Insert Replace	Remov	re Save Cancel

Figure 173. Color Palette Configuration dialog box

- 3. Name the custom color palette.
- 4. Add selected colors to the custom color palette (see "Editing a Custom Color Palette" on page 571).
- 5. Click Save.

The custom color palette appears as the selected palette in the Selected Palette list of the Color Schema view.

6. To apply the new color palette, click **Save Current Settings**.

Deleting Custom Color Palettes

- ✤ To delete a custom color palette
- 1. In the Color Schema view, select the custom color palette in the Selected Palette list.
- 2. Click Delete.

Importing Custom Color Palettes

To import a custom color palette

- 1. In the Color Schema view, click Import.
- 2. Locate the color palette file (XML) and click Open.

The import color palette appears in the Selected Palette list.

Exporting Custom Color Palettes

✤ To export a custom color palette

- 1. In the Color Schema view, select the custom color palette in the Selected Palette list.
- 2. Click Export.
- 3. Name the palette and click **Save**.

Editing a Custom Color Palette

✤ To edit a custom color palette

- 1. Open the custom color palette for editing as follows:
 - a. In the Selected Palette list, select the custom color palette.
 - b. Click Edit.

The Color Palette Configuration dialog box opens.

- 2. Do any of the following:
 - a. To add a color to a custom color palette, select the color in the gradient color chart.
 - b. Click Add.

In the Palette Colors area, the new color appears to the right of the current colors.

-or-

- a. To insert a color in a custom color palette, select its insertion point—the color to the right of the intended position—in the Palette Colors area.
- b. Select the new color in the gradient color chart.

2	olor Palette	Configuration					- • •
Pa	ette name:	Primary Colors					
Pa	ette colors:						
	Advanced o	ptions		Red: Green: Blue: Hue: Saturation: Lightness: Hexcode: Preview:	255 d 0 100% 50% #FF0000		
	Add	Insert	Replace	Remo	ve	Save	Cancel
		Selecte	d color				

Figure 174. Color Palette Configuration dialog box with the basic options view

c. Click Insert.

In the Palette Colors area, the new color appears to the left of the currently selected color.

Olor Palette	Configuration						- • •
Palette name:	Primary Colors						
Palette colors:							
			Red: Green: Blue: Hue: Saturation: Lightness:	255 0 0 0 100% 50%	-		
Advanced o	ptions		Hexcode: Preview:	#FF0000	- 		
Add	Insert	Replace	Remov	/e		Save	Cancel

-or-

- a. To replace a color with another color, select the color to replace in the Palette Colors area.
- b. Select a color in the gradient color chart.

c. Click Replace.

-or-

- a. To remove a color from a custom color palette, select the color in the Palette Colors area.
- b. Click Remove.
- 3. Click Save.

Selecting a Color in the Gradient Color Chart

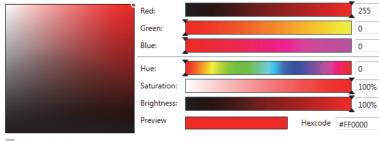
The Color Palette Configuration dialog box that opens when you click either New or Edit in the Color Schema view includes a gradient color chart.

* To select a color in the gradient color chart

Do any of the following:

- Click a color in the hexadecimal color gradient chart.
- Enter an RGB formula or an HSB formula by typing numeric values or using the slider.
- Select the Advanced Options check box and use the separate sliders for red, green, blue, hue, saturation, and brightness.

Figure 175. Advanced Options view



Advanced options

-or-

• Enter the hexadecimal code.

Specifying the Default mzCloud Mass Tolerance Settings

Use the Submit Single Spectrum to mzCloud Options view to set up the mass tolerance settings for a manual mzCloud search.

* To set up the mass tolerance settings for a manual mzCloud search

- 1. In the left pane of the Configuration page, under Miscellaneous settings, select **Submit Single Spectrum to mzCloud Options**.
- 2. Do the following:
 - For the mass tolerance of scans acquired in the Orbitrap mass analyzer, select the units, and then type an appropriate value in the FT Mass Tolerance box.
 - For the mass tolerance of scans acquired in the ion trap mass analyzer, select the units, and then type an appropriate value in the IT Mass Tolerance box.
 - If your processing workflows include the Search mzCloud node, select the **Use** mzCloud Node Settings check box to use the node's settings (see "Search mzCloud Node" on page 209).

3. Click Save the Current Settings.

Table 164 describes the options for submitting single scans to the mzCloud database.

Parameter	Description
FT Mass Tolerance	Specifies the mass tolerance for scans acquired with an FT mass analyzer.
	Default: 12 ppm
IT Mass Tolerance	Specifies the mass tolerance for scans acquired with an ion trap mass analyzer.
	Default 0.4 Da
Use mzCloud Node Settings	When this check box is selected, the application uses the settings in the Search mzCloud processing workflow node if the analysis included this node.
	Default: Selected

Table 164. Submit Single Spectrum to mzCloud Options view

Setting Up a BioCyc Account or Subscription

Follow the instructions in the BioCyc User Login view on the Configuration page to set up your BioCyc subscription or user account. You must have either an organization subscription and a user account or an individual subscription to access all BioCyc databases. A user account alone allows you to access only the EcoCyc database.

To set up a BioCyc account or individual subscription and enter your credentials in the BioCyc User Login view, follow these topics as needed:

- Opening the BioCyc User Login View
- Setting Up a BioCyc Account or Individual Subscription
- Entering, Testing, and Saving Your Account Information

Note To set up a user account or an individual user subscription, you must have Internet access.

Opening the BioCyc User Login View

To open the BioCyc User Login view

1. From the application window, choose **Help > Configuration**.

The Configuration page opens.

- 2. In the left pane, under Miscellaneous settings, select BioCyc User Login.
 - If you have an organization subscription, the Organization Subscription area displays the organization name in green. Otherwise, the area displays the following text in red: No Valid Subscription Found.
 - If you already have a BioCyc user account, entered and tested your credentials, and saved the settings, the Username box displays your email address. Otherwise, the following text appears in red under User Subscription: No Valid Subscription Found.

Figure 176 shows a BioCyc User Login page for a user without a subscription.

Ocompound Discoverer 3.0.0	
<u>File Reporting Lists & Libraries View Window Hel</u>	p
 Server Settings Parallel Options Scratch Directory Options Client Settings Study Management Settings Workflow Editor Settings Color Schema Miscellaneous Settings Submit Single Spectrum to mzCloud Options BioCyc User Login HighChem Fragmentation Options 	BioCyc User Login Credential configuration for BioCyc subscription Organization Subscription Thermo Fisher (organization), Expire: 9/1/2018 5:00:00 AM User Subscription User Subscription User Subscription User Subscription Usersame: Password: Test Credentials No valid subscription found. Subscription Info You must have either an organization subscription and a BioCyc account or an individual subscription to access all BioCyc databases. Access to the EcoCyc database requires only that you create a BioCyc account. To create a BioCyc account, go to https://biocyc.org/preferences.html? status=new. To purchase a BioCyc subscription, go to http:// www.phoenixbioinformatics.org/biocyc/subscriptions.html. You can purchase a subscription for an entire institution or an individual. Once you have created a BioCyc account or have purchased a BioCyc subscription, enter your account credentials in the dialog box shown above. Click Test Credentials to validate your entries. Then, click Save Current Settings to store your account credentials.
	Save Current Settings

Figure 176. BioCyc User Login page for a user without a subscription

Setting Up a BioCyc Account or Individual Subscription

* To set up a BioCyc account or individual subscription

Open the BioCyc User Login view on the Configuration page, and follow the instructions in the view.

Entering, Testing, and Saving Your Account Information

To enter, test, and save your account information

- 1. In the User Subscription area of the BioCyc User Login view on the Configuration page, enter your user name (email address) and password.
- 2. Click Test Credentials.
- 3. Click Save Current Settings.

Figure 177 shows the subscription information for a user who has a user account but does not have an organization subscription or an individual subscription. When running analyses that map compounds to the BioCyc pathways, the user can access only the EcoCyc database.

Figure 177. Settings for a user with a only a user account

(8)	yc User Login ential configuration for BioCyc subscription
	tion Subscription d subscription found.
User Su	scription
Userna	me: bob.chemist@thermofisher.com
Passw	ord: •••••
	Test Credentials Correct password
	Valid credentials: Access to Eco-Cyc
	4

Specifying the Fragmentation Libraries

Use the HighChem Fragmentation Options view to select the fragmentation databases.

Currently, there is only one available fragmentation database. This view is reserved for future use.

* To open the HighChem Fragmentation Options view

In the left pane of the Configuration page, under Miscellaneous Settings, select **HighChem Fragmentation Options**.

14 Setting Up the Configuration Options Specifying the Fragmentation Libraries



Working with the Application Tables

These topics describe the common operations that you can perform on the tables in the Lists and Libraries view, on the pages of a study, and in result files.

Contents

- Moving Up and Down Table Rows
- Sorting Data Tables
- Freezing Table Rows
- Grouping Table Rows
- Changing the Position of Table Columns
- Freezing Table Columns
- Showing or Hiding Table Columns
- Copying Table Entries to the Clipboard
- Filtering the Tables on a Study Page or a List or Library View

Moving Up and Down Table Rows

Use the following procedure to move through the rows in a data table.

* To move up or down through the rows of a result or library table

To move down, press TAB.

To move up, hold down the SHIFT key and press TAB.

Sorting Data Tables

To sort the data tables, see these topics:

- Sorting Table Entries By One or More Columns
- Sorting Table Entries By a Column with a Distribution Map

Sorting Table Entries By One or More Columns

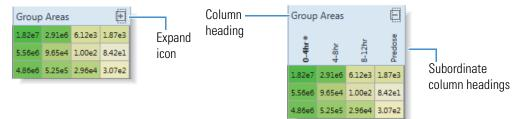
- To sort the rows based on the contents of one or more columns
- 1. Click a column header to sort the rows between ascending order (A, B, C ...) and descending order (Z, Y, X ...), based on the contents of the column.

Note The application treats formulas the same as text strings and sorts them by the order of the characters in the formula string, not by the actual number of elements in the formula.

2. To sort the data by a second column, hold down the CTRL key and click the second column heading.

Sorting Table Entries By a Column with a Distribution Map

- * To sort a table by a column that contains a distribution map
- 1. Click the expand icon to display the vertical headings of the subordinate columns.



2. Select the heading of the subordinate column that you want to sort by.

The selected subordinate column heading appears in bold text.

3. Click the column heading to sort the table rows.

Freezing Table Rows

Use the following procedure to freeze table rows.

✤ To affix rows at the top of the table

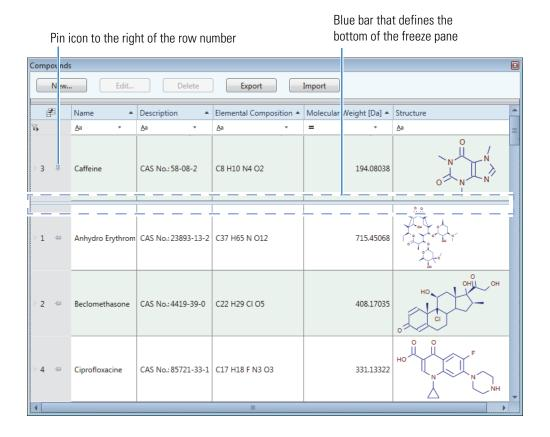
1. Click the pin icon, +, next to the row number of the row that you want to freeze.

The row moves to the top of the table, its pin icon changes to the pinned position, ‡ , and a blue bar that defines the bottom of the freeze pane appears below the fixed row.

2. As you fix additional rows, they move up to the freeze pane in the order selected and their icons change to pinned, [‡]. The row just above the blue bar is the last fixed row.

When you scroll the table, the freeze pane remains at the top. Figure 178 shows a compound library with caffeine in the freeze pane.

Figure 178. Compound library with a freeze pane



Grouping Table Rows

For a table on a study page, use the Enable Row Grouping shortcut menu command to group items by a column heading.

* To group and ungroup the table rows on a study page

1. Right-click the page and choose Enable Row Grouping.

The Group by Area bar appears above the table heading row.

2. Drag the column heading that you want to group by into the Group by Area bar.

Figure 179 shows the column heading for a study factor being dragged to the Group by Area bar.

Stud	ly Definition	Input F	iles Samples Analysis F	Results		
gro	oup by a	area	Drag a field here to group b	y that field	enotype	Column heading
Error	Sample 🔺	File	Sample Identifier	Sample Type	Phenotype	
			I	• •	• •	
	S1	F1	blank	Blank •	n/a •	
	S2	F2	Pooled_ddMS2	Identificatior 🔹	n/a •	
	S3	F3	ZDF_Fatty_1	Sample 🔹	Fatty +	
	S4	F4	ZDF_Fatty_2	Sample *	Fatty •	
	S5	F5	ZDF_Fatty_3	Sample *	Fatty -	
	S6	F6	ZDF_Lean_1	Sample *	Lean -	
	S7	F7	ZDF_Lean_2	Sample 🔹	Lean 🔹	
	S8	F8	ZDF_Lean_3	Sample 🔹	Lean 🔹	

Figure 179. Group by Area bar above the table heading row

3. To ungroup the table rows, drag the column heading out of the Group by Area bar.

Figure 180 shows the study factor (Phenotype) column heading inside the Group by Area bar, and the table rows grouped by the study factor value (Lean, Fatty, or n/a).

Щ	, Add	Files 🛛 👗 F	Remove	File 🔍 Open Contain	ing Folder 🛛 🏶 Ne	ew Analysis 🛛 📢	🗐 Open Analysis Template
St	udy D	efinition	nput Fil	es Samples Analysis	Results		
	F	henotype 🔺	>				
-	Error	Sample 🔺	File	Sample Identifier	Sample Type	Phenotype	
					• •		
9	Fatty	/ (3 items)					
	Error	Sample 🔺	File	Sample Identifier	Sample Type	Phenotype	
		S3	F3	ZDF_Fatty_1	Sample +	Fatty •	
		S4	F4	ZDF_Fatty_2	Sample -	Fatty •	
		S5	F5	ZDF_Fatty_3	Sample +	Fatty +	
9	Lean	(3 items)					
	Error	Sample 🔺	File	Sample Identifier	Sample Type	Phenotype	
		S 6	F6	ZDF_Lean_1	Sample -	Lean 🔹	
		S 7	F7	ZDF_Lean_2	Sample +	Lean •	
		S8	F8	ZDF_Lean_3	Sample +	Lean 🔹	
9	n/a ((2 items)					
	Error	Sample 🔺	File	Sample Identifier	Sample Type	Phenotype	
		S1	F1	blank	Blank *	n/a 🔹	
		S2	F2	Pooled_ddMS2	Identificatior •	n/a •	

Figure 180. Samples table grouped by phenotype

Column heading in the Group by Area bar

4. To turn off the group by row feature, right-click the page and choose **Disable Row Grouping**.

Changing the Position of Table Columns

To change the position of table columns, see the following topics. Changes to tables in any of the Lists & Libraries views are temporary. You can save the layout changes to a result table by applying the File > Save Result View Layout command.

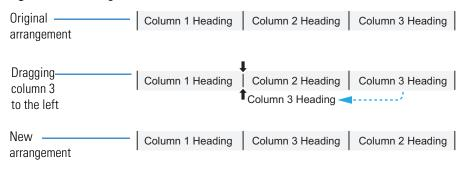
- Changing the Column Order
- Stacking Two Table Columns into One Column

Changing the Column Order

* To change the order of the columns in a library or list table or a result table

To move a column to the left of its current position, drag the column header to the left. Release the mouse button when the cursor () appears over the column delineator (Figure 181).

Figure 181. Moving a column to the left



Stacking Two Table Columns into One Column

To stack two table columns into one column

Drag the column header of the column that you want to stack below the column header of the column that you want on top. Release the mouse button when the cursor ($\rightarrow \leftarrow$) appears over the column heading (Figure 182).

Figure 182. Stacking two columns into one column

Original ———		Column 1 Heading	Column 2 Heading	Column 3 Heading
arrangement	1	Column 1 Entry	Column 2 Entry	Column 3 Entry
	2	Column 1 Entry	Column 2 Entry	Column 3 Entry

Dragging ———		Column 1 Heading		Column 3 Heading
column 3	1	Column 1 Entry	Column 3 Heading Column 2 Entry	Column 3 Entry
into column 2	2	Column 1 Entry	Column 2 Entry	Column 3 Entry

New		Column 1 Heading	Column 2 Heading
arrangement			Column 3 Heading
	1	Column 1 Entry	Column 2 Entry
	1		Column 3 Entry
	2	Column 1 Entry	Column 2 Entry
	2		Column 3 Entry

Freezing Table Columns

In a result table, to more easily compare values in columns that are not next to each other, you can lock columns in place so that they are always visible as you scroll through the unlocked columns.

Note Except for the Input Files table, the Checked column is, by default, the first column in every result table—that is, Checked is the text in first column's heading row.

To lock table columns to the left of the first column

1. Right-click the table and choose Enable Column Fixing.

A pin icon, +, appears to the right of each column heading.

2. Click the pin icons for the columns that you want to move to the left of the first column.

The columns move to the left of the Checked column and their pins face down, Ψ .

Showing or Hiding Table Columns

Use the Field Chooser dialog box to show or hide columns in any of the result tables or tables in the Lists & Libraries views. The changes to tables in any of the Lists & Libraries views are temporary. You can save the layout changes to a result table by choosing Window > Save Layout.

- * To show or hide columns in a library or result table
- 1. Click the **Field Chooser** icon, **P**, in the upper left corner of the table (Figure 183).

Figure 183. Field chooser icon in the upper left corner of the main mzCloud Results table

Field Chooser icon JuzCloud Results ŧ. Checked Structure Name Formula Molecular Weight Best Match 🔻 mzCloud ID KEGG ID 1 -----Caffeine C8 H10 N4 O2 194.08038 97.9 Reference-338 C07481

The Field Chooser dialog box opens with a list of all of the column headers for the current table in alphabetical order (Figure 184).

Field	l Chooser 🛛 🗵
1	Best Match
1	Checked
	Compound Class
1	Formula
1	KEGG ID
1	Molecular Weight
1	mzCloud ID
	mzCloud Library
1	Name
1	Structure

Figure 184. Field Chooser dialog box for the mzCloud Results table

2. In the Field Chooser dialog box, clear the check box for each column that you want to hide. To show those columns again, select their corresponding check boxes.

The table updates and shows or hides your chosen columns immediately.

Copying Table Entries to the Clipboard

You can copy a single table cell, a single table row, or multiple table rows to the Clipboard, and then paste the Clipboard contents into other documents, such as a Notepad text document or Microsoft Office documents.

Note The application does not copy the compound structure in the Structure column of the Expected Compounds library to the Clipboard.

See these topics:

- Copying the Contents of a Single Table Cell to the Clipboard
- Copying the Contents of a Single Row to the Clipboard
- Copying the Contents of Multiple Rows to the Clipboard

Copying the Contents of a Single Table Cell to the Clipboard

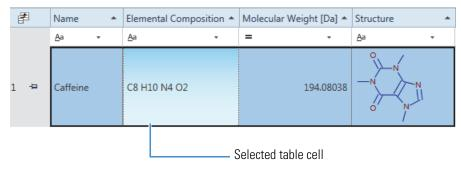
* To copy table cells to the Clipboard

1. Right-click anywhere in the table and choose Cell Selection Mode.

Copy With Headers Copy	Ctrl+C
Clear Selection	
Cell Selection Mode	

2. To select multiple table cells, use the CTRL key or the SHIFT key.

The selected cells turn a lighter blue than the other cells in the row, as shown in the following figure.



3. Right-click and choose **Copy** from the shortcut menu.

Copying the Contents of a Single Row to the Clipboard

To copy a single row to the Clipboard

Do one of the following:

- To copy a single row to the Clipboard, right-click the row and choose **Copy**.
- To copy a single row and the table header, right-click the row and choose **Copy with Headers**.

Copying the Contents of Multiple Rows to the Clipboard

✤ To copy multiple rows to the Clipboard

Do one of the following:

- To copy a range of contiguous rows to the Clipboard, while holding down the SHIFT key, click the first and last row in the range. Then, right-click the last row to open the shortcut menu and choose **Copy** to copy the row contents or choose **Copy** with Headers to copy the row contents and the table header.
- To copy noncontiguous rows to the Clipboard, while holding down the CTRL key, click each row that you want to copy. Then, right-click the last row to open the shortcut menu and choose **Copy** to copy the row contents or choose **Copy with Headers** to copy the row contents and the table header.

Filtering the Tables on a Study Page or a List or Library View

For the tables on the study pages or in the Lists & Libraries views, use the filters in the filter row below the column headers to reduce the number of entries in the current display. The filtering effect is not permanent; closing a filtered list or study removes the filters.

Note For information about filtering the tables in the result files, see "Using Result Filters for Data Reduction and Creating Filter Sets" on page 266.

To set up single-condition filters for the table columns, see these topics:

- Setting Up Single-Condition Filters for the Table Columns
- Setting Up a Single-Condition Wild Card Filter
- Setting Up a Single-Condition Filter for Numeric Data

Related Topics

• Setting Up a Custom Filter with Multiple Conditions

Setting Up Single-Condition Filters for the Table Columns

To set up filters for one or more table columns

- 1. In the table's filter row, do the following for each column that you want to filter by:
 - a. Click the operator symbol (Aa or =) and select an operator from the list.
 - b. Set up the operand by selecting or typing a value in the operand box.

After you set up a filter, the applied filter icon, $T_{\mathbf{x}}$, appears to the right of the operand box, and the table displays only those rows with entries that fulfill the filter condition.

2. To remove a single filter, click the filter icon, $\mathbb{Y}_{\mathbf{x}}$, to the right of the operand box.

Setting Up a Single-Condition Wild Card Filter

To set up a wild card filter for a table column

- 1. In the operator list in the filter row, select Like (Wildcards) or Not Like (Wildcards).
- 2. In the operand box, select or type text and use an asterisk "*" to replace more than one character or use a question mark "?" to replace only one character.

Tip For example, to filter the entries in the transformations library by the presence of nitrogen in the arriving group, do the following in the Arriving Group column:

- Select * Like (Wildcards) in the operator list.
- Type ***N*** in the operand box.

Setting Up a Single-Condition Filter for Numeric Data

* To set up a filter for a table column with numeric data

Do any of the following:

- To set up a filter that uses a specific table entry in the operand list, select any of these operators: Equals, Not Equals, Less Than, Less Than or Equal To, Greater Than, or Greater Than or Equal To.
- To set up a filter that uses any of these operands: (Blanks), (NonBlanks), Above Average, Below Average, Top 10, Top 10 percentile, Bottom 10, or Bottom 10 percentile, select either = Equals or ≠ Not Equals in the operator list.
- To display the top *n* number of entries, select **Top** in the operator list and type an integer value in the operand box.
- To display the bottom *n* number of entries, select **v Bottom** in the operator list and type an integer value in the operand box.
- To display the top *n* percentile of entries, select *** Top Percentile** in the operator list and type a numeric value in the operand box.
- To display the bottom *n* percentile of entries, select **% Bottom Percentile** in the operator list and type a numeric value in the operand box.

Operators and Operands for a Single-Condition Table Filter

The filter for each table column consists of an operator and an operand. In an unfiltered table, the filter row displays the default operator, which is represented by its symbol, and an empty operand box for each column. To set up a table filter, you select the operator from a fixed list, and you select the operand from a list or type a value in the operand box.

The selections in the operator list depend on whether the column contains text or numeric entries. After you select an operator, the operator symbol appears in the filter row to the left of the operand box. For more information about the operator lists, see Table 166 on page 593 and Table 167 on page 594.

For all columns, the operand list includes the following: Custom, Blanks, NonBlanks, and the column entries. For numerical-entry columns, the operand list also includes the following: Above Average, Below Average, Top 10, Top 10 Percentile, Bottom 10, and Bottom 10 Percentile. For more information, see Table 165.

After you set up a column filter, the applied filter icon, $\mathbb{T}_{\mathbf{x}}$, appears to the right of the operand box. Figure 185 shows a filtered Ion Definitions list that reduces the number of displayed entries to 10 by using the total adduct mass. The filter row of the Adducts Total Mass column displays the equals symbol (=) for the mathematical operator, the selection of Top 10 for the operand, and the applied filter icon, $\mathbb{T}_{\mathbf{x}}$.

			Operator				
			Operand				
			Д	pplied	filter icon		
🟦 Lists & Libraries 🗡						~ X	
Ţ	New.	Edit	De ete		Import	Export All	
Expected Compounds		Ion Definition	A dduct s Total Mass	Char	ge	Weight	
85		<u>A</u> a •	= Top 10 -	7 _× =	•	=	F-F
0.0	15 🗁	M+H+ACN	42.033	83	1	50	r
Transformations	2 ⊹⊐	2M+H+ACN	42.033	83	1	50	
	11 🕂	M+2H+ACN	43.041	10	2	50	
28	9 🕂	M-H+FA	44.998	20	-1	50	
Adducts	8 🕂	2M-H+FA	44.998	20	-1	50	
	10 🗁	M-H+HAc	59.013	85	-1	50	
C	9 👳	2M-H+HAc	59.013	85	-1	50	
••	25 🕀	M+Na+ACN	64.015	77	1	50	
Ion Definitions	5 🕀	2M+Na+ACN	64.015	77	1	50	
	16 🕀	M+H+DMSO	79.021	21	1	50	
Da	11 👳	M-H+TFA	112.985	59	-1	50	
Mass Lists							

Figure 185. Ion Definitions list that is filtered by the adduct mass

Table 165 describes the available operand selections and the valid typed operand entries for both text and numeric columns.

Table 165. Operands for the table columns on a study page or Lists & Libraries view (Sheet 1 of 3)

Operand	Description
All table columns	
(Custom)	Applies the custom filter that you set up by using the Custom Filter Selection dialog box.
	A custom filter contains more than one condition. If you set up a single-condition filter, the operand box lists the single condition rather than the (Custom) setting.
(Blanks)	Compatible operators: = Equals and ≠ Not Equals
	= (Blanks)—Displays the table rows that have blank entries in the filtered column.
	≠ (Blanks)—Displays the table rows that have entries in the filtered column.

Operand	Description
(NonBlanks)	Compatible operators: = Equals and ≠ Not Equals
	= (NonBlanks)—Displays the table rows that have entries in the filtered column.
	≠ (NonBlanks)—Displays the table rows that have blank entries in the filtered column.
Selected entry	Table 166 on page 593 describes the compatible operators for text entries. Table 167 on page 594 describes the compatible operators for numeric entries.
	Filters the table rows by using the selected entry and operator.
Typed alphanumeric text or numeric value	Table 166 describes the compatible operators for text entries.Table 167 describes the compatible operators for numeric entries.
	Filters the table rows by using the typed text entry and the selected operator.
Additional selections for	r numeric value columns
Above Average	Compatible operators: = Equals and \neq Not Equals
	= (Above Average)—Displays the table rows with numeric values in the filtered column that are greater than the calculated column average.
	≠ (Above Average)—Displays the table rows with numeric values in the filtered column that are equal to or less than the calculated column average.
Below Average	Compatible operators: = Equals and \neq Not Equals
	= (Below Average)—Displays the table rows with numeric values in the filtered column that are less than the calculated column average.
	≠ (Below Average)—Displays the table rows with numeric values in the filtered column that are equal to or greater than the calculated column average.
Top 10	Compatible operators: = Equals and ≠ Not Equals
	= (Top 10)—Displays the top 10 table rows for the filter condition.
	≠ (Top 10)—Displays the table rows with numeric values in the filtered column that are less than those of the top 10 table rows.

Table 165. Operands for the table columns on a study page or Lists & Libraries view (Sheet 2 of 3)

Operand	Description
Top 10 Percentile	Compatible operators: = Equals and ≠ Not Equals
	= (Top 10 Percentile)—Displays the top 10 th percentile of table rows for the filter condition.
	\neq (Top 10 Percentile)—Displays the table rows with numeric values in the filtered column that are less than those of the top 10 th percentile.
Bottom 10	Compatible operators: = Equals and \neq Not Equals
	= (Bottom 10)—Displays the bottom 10 table rows for the filter condition.
	≠ (Bottom 10)—Displays the table rows with numeric values in the filtered column that are greater than those of the bottom 10 table rows.
Bottom 10 Percentile	Compatible operators: = Equals and ≠ Not Equals
	= (Bottom 10 Percentile)—Displays the bottom 10^{th} percentile of table rows for the filter condition.
	≠ (Bottom 10 Percentile)—Displays the table rows with numeric values in the filtered column that are greater than those of the bottom 10 th percentile.

Table 165. Operands for the table columns on a study page or Lists & Libraries view (Sheet 3 of 3)

Table 166 describes the operators for columns with text entries.

Symbol	Text selection	Effect
=	Equals	Displays the text entries that exactly match the selected or typed operand.
¥	Not equals	Displays the text entries that do not exactly match the selected or typed operand.
<	Less than	For alphabetic text entries, displays the text entries that begin with a letter in the alphabet that comes before the selected or typed operand.
≤	Less than or equal to	_
>	Greater than	-
2	Greater than or equal to	_
	Contains	Displays the text entries that contain the text in the selected or typed operand.
90 1	Does not contain	Displays the text entries that do not contain the text in the selected or typed operand.
*	Like (wildcards)	Displays the text entries that contain the selected or typed text and any additional text represented by an asterisk.
*	Not like (wildcards)	Hides the text entries that contain the selected or typed text and any additional text represented by an asterisk.
H	Match (regular expression)	Displays the text entries that contain the same text as the selected or typed operand.
8-8	Does not match (regular expression)	Displays the text entries that do not contain the same text as the selected or typed operand.
<u>A</u> a	Starts with	Displays the text entries that start with the selected or typed operand.
<u>K</u> a	Does not start with	Displays the text entries that do not start with the selected or typed operand.
a <u>A</u>	Ends with	Displays the text entries that end with the selected or typed operand.
a <u>k</u>	Does not end with	Displays the text entries that do not end with the selected or typed operand.

 Table 166.
 Operators for text columns

Table 167 describes the operators for columns with numeric entries.

Symbol	Text selection	Effect
=	Equals	Displays the numerical entries that equal the selected operand.
¥	Not equals	Displays the numerical entries that are not equal to the selected operand.
<	Less than	Displays the numerical entries that are less than the selected operand.
≤	Less than or equal to	Displays the numerical entries that are less than or equal to the selected operand.
>	Greater than	Displays the numerical entries that are greater than the selected operand.
≥	Greater than or equal to	Displays the numerical entries that are greater than or equal to the selected operand.
*	Тор	Displays the n highest entries in the table, where n equals the integer typed in the operand box.
•	Bottom	Displays the n lowest entries in the table, where n equals the integer typed in the operand box.
%	Top percentile	Displays the entries in the top n^{th} percentile, where n equals the percentage typed in the operand box.
*	Bottom percentile	Displays the entries in the bottom n^{th} percentile, where n equals the percentage typed in the operand box.

Table 167. Operators for numeric columns

Setting Up a Custom Filter with Multiple Conditions

Use the Custom Filter Selection dialog box to set up a custom filter with multiple conditions for a library or list table or a table on a study page.

✤ To set up a custom filter

1. Select (**Custom**) from the operand list for a table column.

The Custom Filter Selection dialog box opens.

Add Condition	- Remove Condition(s)	Group Selected:	🖻 'And' Group	🖺 'Or' Group	Toggle	🖷 Ungroup
Operator		Operand				

Figure 186. Custom Filter Selection dialog box with no conditions

- 2. Do the following for each condition that you want to add to a group:
 - a. Click Add Condition.

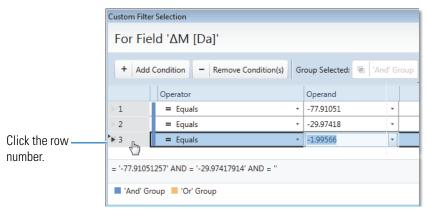
A new table row appears.

b. Select an operator from the Operator list and an operand from the Operand list.

As you add conditions to the group, the application updates the group filter in the gray area below the table.

3. To add the last condition to the group, click its row number.

Figure 187. Clicking the row number in the last row to add the row to the group



The last condition appears in the group filter area. By default, the application applies the AND operator to all of the conditions in the group (Figure 188). A vertical blue bar to the left of the condition rows indicates an AND group.

Custom Filte	er Selection			
For Fie	eld '∆M [Da]'			
+ Add	Condition – Remove Condition	on(s) Gr	oup Selected:	🖻 'And' Group
	Operator		Operand	
▶ 1	= Equals	*	-77.91051	•
≥ 2	= Equals	•	-29.97418	•
₩ 3	= Equals	-	-1.99566	.
	1257' AND = '-29.97417914' AND = roup 📕 'Or' Group	-1.99566	357'	
				- Third condit

Figure 188. Group filter with three conditions and the AND group operator

4. To change the group operator from AND to OR or from OR to AND, click Toggle.

An orange bar to the left of the condition rows indicates an OR group.

Figure 189. Group filter with three conditions and the OR group operator

	Operand
•	-77.91051 🔹
•	-29.97418 *
-	-1.99566 •

- 5. To add an overlapping group to the filter, do the following:
 - a. Select the rows that you want to group, using the SHIFT key for contiguous rows or the CTRL key for noncontiguous rows.

The selected rows are highlighted in blue and the 'And' Group and 'Or' Group buttons become available.

Figure 190.	Selection	of two	noncontiguous rows

+ Add	Condition - Remove Con	dition(s) Group Selected: 🖷 'An	nd' Group 🛛 🕒 'Or' Group
	Operator	Operand	
1	= Equals	-77.91051	•
2	= Equals	-29.97418	•
3	= Equals	-1.99566	.
€ 4	= Equals	· 42.01056	*

b. Specify the group type by clicking 'And' Group or 'Or' Group.

The application applies the second group definition and the Ungroup button becomes available.

Figure 191. A set of filter conditions with two groups

+ Add	Condition Remove Condition(s) G	roup Selected: 🖭 'And' Group	🖫 Or' Group	🔀 Toggle 🗮 Ungroup
	Operator	Operand		
1	= Equals	-77.91051	•	
2	= Equals	-29.97418	•	
3	= Equals	42.01056	-	
> 4	= Equals	-1.99566	•	

- 6. To remove conditions from a group, select the conditions and click Ungroup.
- 7. To apply the filter, click **OK**.

The Custom Filter Selection dialog box closes, the text (Custom) appears in the operand box, and the application applies the custom filter to the entries in the selected filter column.

Custom Filter Selection Dialog Box

Table 168 describes the features of the Custom Filter Selection dialog box.

Table 168. Custom Filter Selection dialog box features

Feature	Description
Buttons or icons	
+ Add Condition	Adds a blank condition row to the condition table.
- Remove Condition(s)	Removes the selected conditions. Selected conditions are highlighted in blue.
'And' Group	When the filter contains more than one group, applies the AND group type to a set of selected conditions.
'Or' Group	When the filter contains more than one group, applies the OR group type to a set of selected conditions.
Toggle	Changes the selected AND group to an OR group or the reverse.
Ungroup	As you create groups, group label columns appear to the left of the Operator column.
	When conditions belong to more than one group, removes the second group condition for the selected conditions.
ОК	Closes the dialog box and applies the filter conditions.
Cancel	Closes the dialog box without applying the filter conditions.
Table	
Operator column	Use to select an operator for the filter condition.
	See Table 166 on page 593 for a list of the operators for the text entry columns. See Table 167 on page 594 for a list of the operators for the numerical entry columns.
Operand column	Use to select or type an operand for the filter condition.
	See Table 165 on page 590 for a list of the operands for the library columns.
Third column	Displays comments about the filter condition. For example, this box displays "Condition is empty" until you define the operator and the operand for a condition.
Filter description area	
This area, which is highlig	hted in gray, displays the group filter conditions.



Working with Molecular Networking

Use the new molecular networking (beta) feature to investigate possible relations between compounds in the Compounds table.

Contents

- Using the Molecular Networking Feature
- How the Generate Molecular Networks Node Works
- Information Displayed in the Similar Compounds Table
- Exporting the Molecular Network to the Viewer
- Working with the Molecular Network Viewer

Using the Molecular Networking Feature

To use the molecular networking feature, follow this process:

- 1. Process your raw data files with a processing workflow that includes the Generate Molecular Networks node, which is a node under Compound Scoring.
- 2. To review the processed results, open the result file.
- 3. (Optional) Using the Result Filters view, apply a filter to display only the compounds of interest.
- 4. Do one or both of the following:
 - To view a table of similar compounds, in the Compounds table, select a compound of interest. Then, open its related Similar Compounds table.
 - -or-
 - Export the molecular network to the viewer.

Related Topics

- How the Generate Molecular Networks Node Works
- Information Displayed in the Similar Compounds Table

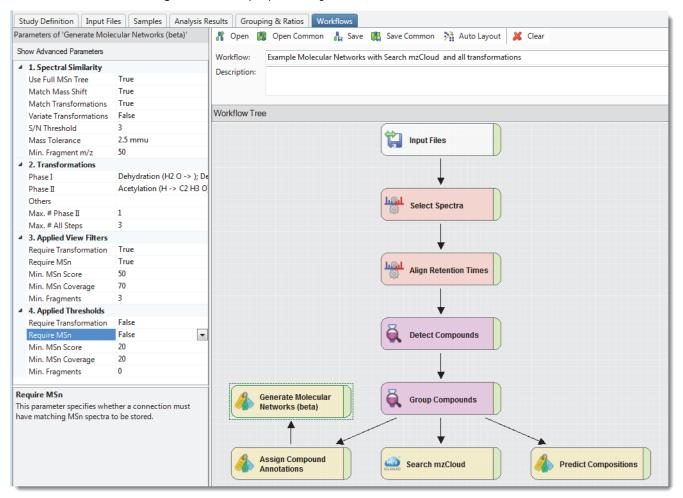
- Exporting the Molecular Network to the Viewer
- Using Result Filters for Data Reduction and Creating Filter Sets

How the Generate Molecular Networks Node Works

To express the similarity of a pair of compounds, the Generate Molecular Networks node uses the assigned elemental compositions and fragmentation data. The composition information is used to reject pairs having unexpected elemental composition differences, while the best MSn fragmentation scans for both compounds are compared by using the FISh scoring algorithm to indicate their structural similarity.

A minimal processing workflow that includes the Generate Molecular Networks node consists of untargeted compound detection, grouping, and annotation assignment. To validate the elemental composition differences, the workflow must contain at least one search node or the Predict Compositions node.

Figure 192. Example processing workflow with the Generate Molecular Networks node



The input to the Generate Molecular Networks node is a list of annotated compounds, and the output from the node is a table of similar compounds for each detected compound.

Note If you do not select any transformations for the Generate Molecular Networks node, the node does not consider the elemental compositions of the detected compounds for scoring.

The node processes the data as follows:

- 1. Generates possible transformation pathways according to the user-specified settings.
 - If there are multiple pathways to the same elemental composition, the node selects the shorter pathway.
 - If there are compounds with the same elemental composition (isomers with different retention times), the node adds a pathway of length 0.
- 2. For each pair of compounds that have assigned elemental compositions, the node finds a matching pathway that explains the elemental composition difference between the two compounds.
- 3. For each pair of compounds that have fragmentation data, the node calculates the spectral similarity by FISh Scoring as follows:
 - a. Matches the fragments for both compounds by directly comparing their masses.
 - b. For the remaining unmatched fragments, the node does the following to match the fragments:
 - Uses the mass shift of the assigned pathway to match "shifted" fragments (if enabled). In addition, it uses all possible permutations of the individual pathway steps (if enabled).
 - Uses the mass shift between the two compounds to match "shifted" fragments (if enabled).
 - c. Calculates the similarity scores.
- 4. Applies specified rules and thresholds to the connections (matched pairs).
- 5. Stores valid connections to the results file.
- 6. Applies specified view filters on the results table.

Related Topics

• Similar Compounds (Beta) Related Table

Information Displayed in the Similar Compounds Table

The Generate Molecular Networks node stores all the valid connections between each pair of compounds and creates a related table of Similar Compounds. The table shows all the connections between the selected compound and its related compounds, with additional information about the similarity between them.

You can consider each stored connection as a reaction, where one compound is a substrate that is converted into a product through a specific transformation pathway. The Direction column indicates the direction of the reaction between the selected compound in the Compounds table and the similar compound in the Similar Compounds table.

A Forward connection indicates that the selected compound is a substrate, to which the transformation has been applied to generate the similar compound. A Reverse connection indicates that the selected compound is a product of applying the transformation to the similar compound.

In Figure 193, paraxanthine is listed as a demethylation product of caffeine (substrate).

Figure 193. Caffeine and one of its demethylation products

•	Compo	unds C	Compounds	per File Feat	ures mzCloud Re	ults Input Files										
	1	Checked	Name		Formula	Annotation	Source 🛨	FISh Coverage	Molecular V	Veight	RT [min]	Area (Max.) 🔻	# mzCloud Results	mzCloud Best Mate	ch MS2	Area 🛨
2	4		Caffeine		C8 H10 N4 O2				194.0	08026	3.780	3425685171	3	95.	4	3.43e9
3	4		5-Hydroxy	romeprazole	C17 H19 N3 O4 S				361.1	10917	4.919	1781181583	2	92.	9	1.78e9
Hide Related Tables																
Str	ucture F	Proposals	Compo	unds per File	Predicted Compositi	ons Similar Compo	ounds (beta)	😵 mzCle	oud Results							
	ŧ.	Checked	Direction	Mass Shift [Da]	Composition Chan	e Transformations	Transforma	ation Mass [Da]	# Fragments	s MSn S	Score For	ward Cov. [%]	# Forward Matches	Reverse Cov. [%] #	‡ Revers	e Matches
1	4		Forward	-14.01565	C-1 H-2	demethylation		-14.01565	6	i -	67	67	10	67		4
2	. ÷		Forward	-14.01565	C-1 H-2	demethylation		-14.01565	6		64	62	8	67		4
3	÷ +=		Forward	0.00000		Isomer		0.00000	6		78	89	8	67		4
Hide Related Tables Compounds																
	ŧ.	Checked	Direction	Name	Formula	Annotation Source 🗄	FISh Cove	erage Molecul	ar Weight RT	[min]	Area (Ma	k.) ▼ # Fragme	ents # mzCloud Res	ults mzCloud Best	Match	MS2 Area 🛨
1	-12		Forward	Caffeine	C8 H10 N4 O2			1	94.08026	3.780	3425685	5171	6	3	95.4	3.43e9
2			Reverse	Paraxanthine	C7 H8 N4 O2			1	80.06455	3.214	1066578	3149	15	6	93.7	1.07e9

Exporting the Molecular Network to the Viewer

After you process a set of input files with a processing workflow that includes the Generate Molecular Network node, you can export the molecular network for a specified number of compounds to the molecular networking viewer.

✤ To export the molecular network to the viewer

- 1. (Optional) To display only the compounds of interest, filter the Compounds table.
- 2. (Optional) Sort the Compounds table by the columns of interest.

Note The application only exports compounds that appear in the table in the order that they appear in the table, beginning with row 1. It does not export compounds that are hidden by applied result filters or compounds in row numbers greater than the Compounds Limit value.

3. Right-click the Compounds table and choose Molecular Networks (Beta).

The Export Molecular Networks (Beta) dialog box opens with the following default settings:

- Destination: Folder where the result file resides
- Name: Name of the result file
- Compounds Limit: 500
- Open Viewer After Export: Clear
- 4. Do any of the following:
 - To change the destination folder, click the browse icon and select another folder.
 - To change the name of the final folder where the application stores the web page components, type a different name in the Name box.
 - To change the maximum number of compounds to export, type a number from **1** to **1000** in the Compounds Limit box.

Note Increasing the number of compounds increases the processing time.

- To automatically open the viewer in the default browser after clicking Export, select the **Open Viewer After Export** check box.
- 5. Click Export.
- 6. If you did not select the Open Viewer After Export check box, open the named destination folder. Then, right-click the index.html file and choose the web browser of interest.

The viewer opens in the selected or default web browser.

Working with the Molecular Network Viewer

Use the molecular network viewer to visualize the similarity between the compounds of interest in the Compounds table.

When you export compounds to the molecular networking viewer, the application processes the data and displays a molecular network simulation with the filter and threshold settings that you specified in the Generate Molecular Networks workflow node.

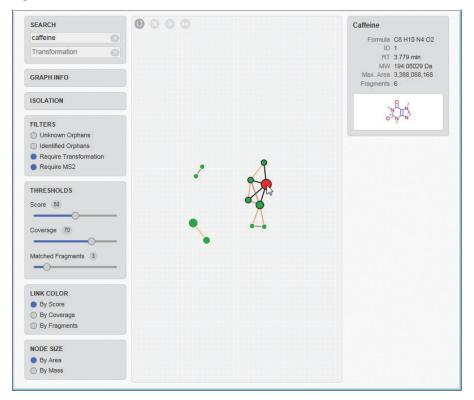
In the left pane of the viewer, you can change the settings for the filters and the thresholds, change the basis of the link color and the node size, search for a named compound or transformation, and isolate the display to a specific cluster.

In the right pane of the viewer, you can view information for each node or link that you point to in the graph area.

See these topics:

- Color-Coding for the Nodes
- Sizing a Node By the Compound's Peak Area or MW
- Colorizing a Link By Its Score, Coverage, or Number of Fragments
- Interactive Functions in the Molecular Network Viewer
- Search Pane

Figure 194. Molecular network viewer with a cluster for caffeine



Color-Coding for the Nodes

In the molecular network viewer, the color of a compound's node indicates the identification confidence.

Table 169. Node colors

Color	Meaning
Light green	The workflow identified the compound by its fragmentation spectra during an mzCloud or mzVault library search.
Dark green	The workflow identified the compound by its formula and a database match.
Blue	The workflow only determined the compound's formula.
Gray	The workflow only determined the compound's mass.

Sizing a Node By the Compound's Peak Area or MW

The node size is proportional to the Max. Area or MW of a compound, depending on the option you select.

- To select whether the node size is proportional to the compound's peak area or MW
- 1. In the left pane of the molecular network viewer, open the Node Size area.
- 2. Select one of these options: By Area or By Mass.

Colorizing a Link By Its Score, Coverage, or Number of Fragments

The color of a link is displayed in the range of gray > orange > dark-red and is proportional to the MSn Score, the Forward or Reverse Coverage (max), or the number of Forward or Reverse Matches (max), depending on the option you select. If a connection has no fragmentation data, the link appears as a gray-dashed line.

✤ To select a different option to colorize a link by

- 1. In the left pane of the molecular network viewer, open the Link Color area.
- 2. Select one of these options: By Score, By Coverage, or By Fragments.

Tip By default, the link color is a function of the MSn score. To investigate the clusters, try colorizing the links by the other options.

Note The length of each link has no meaning.

Interactive Functions in the Molecular Network Viewer

Table 170 describes the interactive tasks that you can perform in the molecular network viewer.

Table 170. Interactive functions in the molecular network viewer
--

Task	Action
Show information about a compound.	Point to the compound.
(graph node)	Information about the compound appears in the right pane, and all the compound's relations are highlighted in the graph area.
Show information about a graph link	Point to the link (connecting line).
(relation).	Information about the connection between the two compounds appears.
Zoom in or out of an area of the graph.	Point to the area of the graph where you want to change the zoom level, and then use the mouse wheel.
There are four icons in the upper left of	the graph.
Reset the pan and zoom.	Click the Reset Pan and Zoom icon, 🕣 . Or, double-click anywhere in the graph.
Move a cluster.	Drag one of its nodes.
Refresh the simulation.	Click the Refresh icon, 🔘 .
Stop the simulation process.	Click the Stop Simulation icon, 🗙 .
Enter or exit the isolation mode.	Double-click a node to isolate the node and its connections, and then click the Isolation mode icon, 😒, to exit the isolation mode.

Search Pane

Use the Search pane of the molecular network viewer to highlight nodes or links in the graph area by searching for the name of a compound or the name of a transformation, respectively.

To do the following	Do this
Highlight nodes, in red, for compounds with a specified text string in their names.	In the Search pane, type the compound's name in the Compound box. The viewer highlights all compounds with the
	specified text string in the graph area.
Highlight links, in red, for the same named transformation.	In the Search pane, type the transformation's name in the Transformation box.

Graph Info Pane

The Graph Info pane of the molecular network viewer provides information about the exported data. Each bar shows relative portion of currently visible items in blue (for example, the number of visible links with fragmentation data) as well as relative portion of all items of particular type in dark gray (for example, the number of all links with fragmentation data). Finally, the light-gray background of a bar represents the total number of items (for example, the number of all the links). The tooltip provides the exact numbers if needed.

Parameter	Description
Nodes	Indicates the number of nodes in the graph area.
Orphans	Indicates the number of nodes without any connection (according to the current filters and thresholds).
Identified	Indicates the number of compounds with an assigned name.
Unknown	Indicates the number of compounds without an assigned name.
Links	Indicates the number of all connections.
Transform.	Indicates the number of connections with assigned transformations.
MS2	Indicates the number of connections with fragmentation data.
Both	Indicates the number of connections with assigned transformations and fragmentation data.

Table 171. Graph Info pane of the molecular network viewer

Isolation Pane

If the main graph is in the Isolation mode, the Isolation panel provides dynamic control over the maximum allowed depth of the graph (see below for more details).

Filters Pane

Use the filters in the Filters pane of the molecular network viewer to limit the amount of visible data.

Table 172. Filters pane (Sheet 1 of 2)

Filter	Description
Unknown Orphans	Select to display compounds without a name or connection.
	Default: Not enabled
Identified Orphans	Select to display named compounds without a connection.
	Default: Not enabled

Table 172. Filters pane (Sheet 2 of 2)

Filter	Description
Require Transformation	Clear to display links without assigned transformations.
	Default: Enabled
Require MS2	Clear to display links without fragmentation data.
	Default: Enabled

Thresholds Pane

To show or hide low confidence relationships between compounds, adjust the settings in the Thresholds pane. By default, the thresholds are set to those specified in the processing workflow.

Threshold	Description			
Score	Specifies the minimum MSn Score for a connection to be visible.			
	Range: 0 to 100			
Coverage	Specifies the minimum Forward or Reverse Coverage for a connection to be visible.			
	Range: 0 to 100			
Matched Fragments	Specifies the minimum number of matched fragments for a Forward or Reverse search for a connection to be visible.			
	Range: 0 to 20			

Compound or Transformation Information

Each time you point to a link or node, available information appears on the right side of the viewer. For a link, the information pane first displays the assigned pathway name, elemental composition, mass difference and fragmentation scores, followed by the reaction from one compound to the other. For each compound, it shows the assigned name, elemental composition, retention time, molecular weight, max. area, and number of used fragments.

The information automatically disappears from the right pane when you move the pointer away from the node or link.

Tip To prevent the information about the selected node or link from disappearing or changing when you move the mouse pointer, hold the SHIFT key while you move the pointer.

Using the Isolation mode

Use the isolation mode to display only the cluster of interest. The graph displays the connections to the selected node up to the maximum reaction depth that you specify under Visible Depth.

* To turn on the isolation mode to isolate a specific cluster

- 1. Double-click the node of interest.
- 2. To exit the isolation mode, click the **Toggle Isolation** icon, S, or double-click the orange node—that is, double-click the node that you clicked to enter the isolation mode.

16 Working with Molecular Networking Working with the Molecular Network Viewer

17 -

Testing Communication to the Online Databases

A typical Compound Discoverer analysis that identifies unknown compounds searches mass spectrum databases on the Internet. To run these searches, the application must have unblocked access to the mass spectral databases on the Internet.

Follow these topics to test and troubleshoot the application's access to the online mass spectrum databases.

Contents

- Running the Communication Tests
- Checking the URLs for the Online Databases in Your Browser
- Specifying the IP Address of the Proxy Server
- Setting the Correct Time and Time Zone on the Processing Computer

To test and troubleshoot the application's access to the online databases

1. Run the communication tests (see "Running the Communication Tests" on page 612).

If the communication tests succeed, the application has access to the online databases.

- 2. If a communication test fails, do the following as needed:
 - If only the mzCloud communication test fails, check the Date and Time settings on the processing computer (see "Setting the Correct Time and Time Zone on the Processing Computer" on page 615).
 - If the Check Subscription test for the BioCyc database fails, check the subscription information in the BioCyc User Login view of the Configuration page.

IMPORTANT If you do not have an organization subscription for the BioCyc database, you must create a BioCyc user account or obtain and individual subscription, and then enter, test, and save your account credentials in the BioCyc User Login view (see "Setting Up a BioCyc Account or Subscription" on page 575).

• If any of the other communication tests also fail, check the access to the URLs for the online databases (see "Checking the URLs for the Online Databases in Your Browser" on page 613).

If you can access the URLs for the online databases through your browser, but the communication tests still fail, the firewall or proxy setting for your company network is blocking the application's access to the online databases.

- 3. If the communication tests fail, but you can access the URLs for the online databases, do the following as needed:
 - If a firewall is blocking the application's access to the online databases, ask your IT department to make sure that the company firewall is not blocking "Compound Discoverer" or "Compound Discoverer Server" from accessing the URLs. The application uses the following protocol: http port 80 and .
 - If a proxy setting is blocking access, see "Specifying the IP Address of the Proxy Server" on page 614.

Running the Communication Tests

Use the Communication Tests dialog box to test your processing computer's access to the online databases.

- * To verify that your computer has access to the external databases
- 1. From the menu bar, choose **Help > Communication Tests**.

The Communication Tests dialog box opens.

- 2. To open the page for the database that you want to access, click its tab.
- 3. Click Run Tests.

Figure 195 shows the communication tests in progress.

Figure 195. BioCyc co	ommunication tests
-----------------------	--------------------

🖉 Communication Tests	x
BioCyc KEGGPathway mzCloud ChemSpider	
Checking service availability Succeeded	
Succeeded	
Check Subscription Succeeded	
Checking formula search Succeeded	
Checking MW search Succeeded	
BioCyc: Checking formula search - Performing test search by formula 'C6H6'	*
-> Succeeded	
Succeeded	4 III
Clear Results Copy Info Run Tests Close	

4. If the tests are successful, your computer has access to the required databases on the Internet. If only the mzCloud test fails, check the Date and Time settings for the processing computer ("Setting the Correct Time and Time Zone on the Processing Computer" on page 615). If any of the other tests also fail, check the access to the URLs in your browser (see "Checking the URLs for the Online Databases in Your Browser" on page 613).

Checking the URLs for the Online Databases in Your Browser

Table 174 lists the URLs for the online mass spectrum databases. If a communication test fails, test the URL for the affected database.

Database	URL
mzCloud Identity	https://identity.mzcloud.org/
	https://www.mzcloud.org/Services/MzCloudApiV1.svc
	https://www.mzcloud.org/Services/MzCloudApiLightService.svc
ChemSpider	http://www.chemspider.com
	http://chemspider.com/MassSpecAPI.asmx
	http://chemspider.com/Search.asmx

Table 174. URLs of online mass spectrum databases (Sheet 1 of 2)

Database	URL
KEGG: Kyoto	http://www.kegg.jp/
Encyclopedia of Genes and Genomes	http://rest.kegg.jp
BioCyc	https://biocyc.org/
	https://biocyc.org/web-services.shtml

Table 174. URLs of online mass spectrum databases (Sheet 2 of 2)

Specifying the IP Address of the Proxy Server

If the communication tests fail but you can access the online databases through your browser, follow this procedure to specify the IP address of the proxy server.

* To configure the IP address of the proxy server

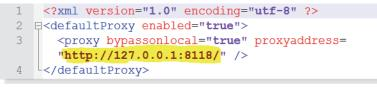
- 1. Go to *drive*:\Program Files\Thermo\Compound Discoverer 3.0\bin\Config.
- 2. Open the **Proxy.config** file in Notepad.
- 3. Remove the text that is highlighted in yellow in Figure 196—that is, remove the XML comment delimiters: <!-- and -->.

Figure 196. Proxy configuration setting with XML comment delimiters

1	xml version="1.0" encoding="utf-8" ?
2	<pre> G<defaultproxy enabled="true"> </defaultproxy></pre>
3	· ¢ </th
4	<pre><pre><pre><pre><pre>proxy bypassonlocal="true" proxyaddress="http://127.0.0.1:8118/" /></pre></pre></pre></pre></pre>
5	>
6	

4. Replace the text that is highlighted in yellow in Figure 197 with your company proxy address.

Figure 197. Default proxy address highlighted in yellow



Setting the Correct Time and Time Zone on the Processing Computer

The mzCloud communication test includes a validation of the date and time settings on the processing computer. If the mzCloud communication test fails, but the other communication tests succeed, check the date and time settings for the processing computer.

- * To check the time and time zone settings
- Open the Date and Time dialog box. For the Windows 7 operating system, open the Control Panel. In the View By list, select Category, and then choose Clock, Language, and Region > Date and Time.

💼 Date and Time
Date and Time Additional Clocks
Date: Saturday, June 03, 2017 Time: 4:35:58 PM
(UTC-08:00) Pacific Time (US & Canada) Change time <u>z</u> one
Daylight Saving Time ends on Sunday, November 05, 2017 at 2:00 AM. The clock is set to go back 1 hour at that time.
☑ Notify me when the clock changes
Get more time zone information online How do I set the clock and time zone?
OK Cancel Apply

2. Make sure that both the date and time and time zone settings are correct.

Tip If your computer is not part of a network domain that synchronizes the computer's clock to the network server, you can use an Internet server to synchronize the computer's clock.

3. If the Internet Time tab is available, click it and synchronize the computer's clock with an Internet server.

17 Testing Communication to the Online Databases Setting the Correct Time and Time Zone on the Processing Computer

Experiment Design for Comparison Statistics

To understand how and when to use the biological replicate study factor, see these topics.

Contents

- Biological Versus Technical Replicates
- Non-Nested Versus Nested Experiment Designs

Biological Versus Technical Replicates

Biological replicates are samples from biological individuals (or non-biological entities) of the same type under the same conditions and provide a measure of the variability associated with these conditions.

Technical replicates are replicate samples from the same entity under the same conditions. Technical replicates from the same entity under the same conditions provide a measure of the sampling error, and replicate injections from the same sample solution provide a measure of the instrument error.

You can add only one biological replicate factor to a study. The application treats study factors nested under a biological replicate factor as technical replicates.

Non-Nested Versus Nested Experiment Designs

When you add a biological replicate factor to study, you can set up two different experiment designs—nested and non-nested. In non-nested experiments, the biological replicates are independent of each other—that is, you do not reuse individual entities to study multiple condition states (study factor items).

In non-nested experiments, different sets of biological replicates are used for each condition. Figure 198 shows a non-nested experiment where replicate samples are taken from two sets of rats under two conditions. Tom, Jerry, and Hector are fed a normal diet; and Tai, Fernando, and Pierre are fed a low-calorie diet.

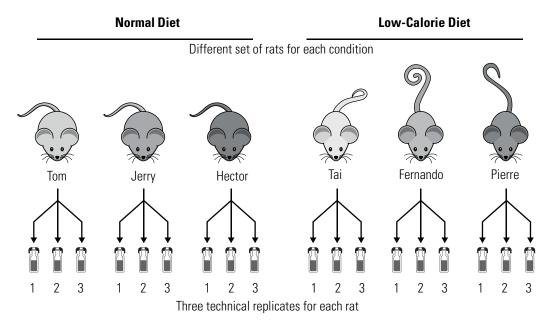


Figure 198. Non-nested design with independent sets of rats for the two dietary conditions

In nested experiments, the same set of biological replicates are used for each condition. Figure 199 shows a nested experiment where replicate samples are taken from the same three rats under two conditions—a normal diet and a low-calorie diet. Tom, Jerry, and Pierre are the biological replicates.

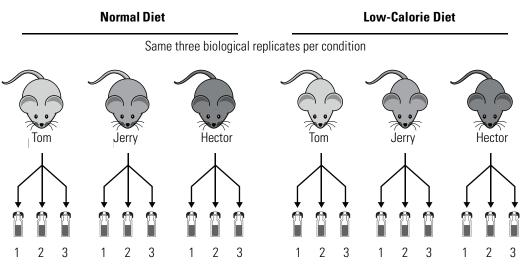


Figure 199. Nested design with the same three rats under two dietary conditions

Three technical replicates for each rat

To set up the comparison ratios for a nested design with technical replicates, add the following study factors:

- A factor for the variable being studied with an itemized list of the variable states
- A biological replicate factor with an itemized list of the entities being studied
- (Optional) A factor for the technical replicates

Figure 200. Study factors for the nested design experiment

Figure 200 shows the study factors for the experiment shown in Figure 199.

Paste Copy Add -Study Factors Numerical factor Injection replicate Edit × 1 2 3 Categorical factor B Diet Edit X Low-cal Diet Normal Diet **Biological Replicate factor** 🎎 Rat Edit × Jerry Pierre Tom

Figure 201 shows the generated sample groups and ratios for the nested design. The sample groups—Low-cal Diet and Normal Diet—are highlighted in green. The two sample groups contain the same values for the Rat Biological Replicate factor—Tom, Jerry, and Pierre. The technical replicates for each biological replicate are grouped together and the biological replicates are highlighted in yellow. As shown in the Generated Ratios area, for each selected denominator, the Differential Analysis node calculates one group ratio and individual ratios for each biological replicate. In the result file, the Compounds table includes a Ratio column for the group ratio and Bio. Rep. Ratio columns for the biological replicate ratios.

Note The application calculates p-values as follows:

- Uses the t-test when comparing two sample groups.
- Uses ANOVA when comparing more than two sample groups.

mple Group and Ratio Specification	Generated S	ample Groups
Study Variables	Low-cal D	iet
🔲 File		Sample Jerry Low-cal Diet 1 F1: Low-cal Diet_Jerry_Rep 1
Diet	Jerry	Sample Jerry Low-cal Diet 2 F2: Low-cal Diet_Jerry_Rep 2
Injection replicate		Sample Jerry Low-cal Diet 3 F3: Low-cal Diet_Jerry_Rep 3
_		Sample Pierre Low-cal Diet 1 F4: Low-cal Diet_Pierre_Rep 1
Sample Type	Pierre	Sample Pierre Low-cal Diet 2 F5: Low-cal Diet_Pierre_Rep 2
Variables printed in italics contain only a single value.		Sample Pierre Low-cal Diet 3 F6: Low-cal Diet_Pierre_Rep 3
		Sample Tom Low-cal Diet 1 F7: Low-cal Diet_Tom_Rep 1
Manual Ratio Generation	Tom	Sample Tom Low-cal Diet 2 F8: Low-cal Diet_Tom_Rep 2
Numerator:		Sample Tom Low-cal Diet 3 F9: Low-cal Diet_Tom_Rep 3
Numerator: Add Ratio	Normal Di	et
Denominator:		Sample Jerry Normal Diet 1 F10: Normal Diet_Jerry_Rep 1
	Jerry	Sample Jerry Normal Diet 2 F11: Normal Diet Jerry_Rep 2
Bulk Ratio Generation		Sample Jerry Normal Diet 3 F12: Normal Diet_Jerry_Rep 3
Denominators to be used:		Sample Pierre Normal Diet 1 F13: Normal Diet_Pierre_Rep 1
Diet : Low-cal Diet Image: Operating the second s	Pierre	Sample Pierre Normal Diet 2 F14: Normal Diet_Pierre_Rep 2
		Sample Pierre Normal Diet 3 F15: Normal Diet_Pierre_Rep 3
		Sample Tom Normal Diet 1 F16: Normal Diet_Tom_Rep 1
	Tom	Sample Tom Normal Diet 2 F17: Normal Diet_Tom_Rep 2
		Sample Tom Normal Diet 3 F18: Normal Diet_Tom_Rep 3
	Generated R	atios (Nested Design) 🛛 💥 Clear
	X Low-ca	I Diet / Normal Diet
	Low-	cal Diet Jerry / Normal Diet Jerry
	Low-	cal Diet Pierre / Normal Diet Pierre
Add Ratios		cal Diet Tom / Normal Diet Tom

Figure 201. Nested design example

Index

A

Abort command, Job Queue 128 Activate button, License Manager 554 activation code, licensing 555 activation request (.req) file 558 activation response file (activation.xml) 557 Activation Type parameter 171, 179 Actual Size icon, reporting 468 Add Column command 523 Add Files command, study 101 Add Items icon, reporting 467 Add Plot command 302 Add Structure Proposal command 279 adding adducts to the library 503 biological replicate factors to a study 68, 81 categorical factors to a study 79 charges to a structure 540 column to a mass list 523 compounds to the library 501 custom explanations to a result file 260 fragments lists to the Compound Classes list 30 input files to a new study 75 input files to the Analysis pane 118 ion definitions to the library 508 manual peaks to a specialized trace 299 mass lists to the library 518 numeric factors to a study 80 page break after the page footer of a report 462 transformations to the library 512 Adduct Editor 503 adducts deleting 503 editing 504 library 502 table, hidden 257 Align Columns icon, reporting 467

Align Retention Times node limitations 174 using 166 Alignment Model parameter 166 Alignment property, reporting 475 alignment, chromatographic 18, 25, 174, 365 Allow Aromatic Cleavage parameter 265 ammonium adduct 192 analyses reprocessing 131 reviewing 130 sample groups and ratios for 92, 120 starting 60 submitting to the job queue 122 troubleshooting 57, 120 Analysis pane 60, 133 Analysis Results page of a study 129 Analysis Sequence Details window 130 analysis summary 275 Analysis Validation Issue prompt 60 Analysis Validation Issues prompt 122 Analyze Labeled Compounds node 189 AnchorBottom property, reporting 475 AND logical conjunction, filtering with 271-272, 274 Angle property, reporting 478 Annotate Full Spectrum Tree check box 265 annotations copying 310 editing 259 fragmentation scan 307 Appearance properties, reporting 474 Appendix section, reporting 466 application menu bar 7 toolbar 12 applications, installing Compound Discoverer xxii mzVault xxiv

Index: B

Apply Dealkylation parameter 237 Apply Dearylation parameter 237 Apply Filters button, Result Filters view 274 Apply FISh Scoring check box 261 Apply mzLogic node 223 arranging windows, tool for 45 Arriving Group data entry box 513 parameter 514 result table column 400 Transformations library column 512 arrows layout guide tool 45 Metabolika pathway 531 As Batch check box, Analysis pane 122, 133 Assign command, Input File Characterization page 84 Assign Compound Annotations node 200 Atom Properties dialog box using for charged fragments 540 using for neutral compounds 551 Auto Layout command, Workflows page 138 AutoReplaceFields property, reporting 478 auto-save option for studies 567 Autosize property, reporting 478 Average Peak Width parameter 19, 232 Avg. Exchange column 407 axis labels Hierarchical Cluster Analysis view 375 Histogram Chart 317 Scatter Chart 328, 330

B

BackColor property, reporting 475 Bar Chart page, displaying 319 Barcode design item 472 BarHeight property, reporting 475 base ions 29, 192 batch normalization 155 Behavior properties, reporting 478 best fit MS scan 394 best MS1 scan for a compound 33 best MS2 scan for a compound 33 BioCyc credentials, saving 576 BioCyc Pathways table 425 BioCyc Pathways view 362 BioCyc Results table 425, 428 BioCyc User Login view 575 Biological Replicate Factor 68, 81, 104, 617 blanks Blank check box for sample type 254

Blank sample type 69, 84, 112 Exclude Blanks parameter 199 Mark Background Compounds node 21, 26, 240 Normalize Areas node 199 blocked IP addresses 611 blue circles, KEGG database 360 bond multiplicity, changing 550 borders, adding to design items 470 bracket, using in elemental composition formula 144 Bulk Ratio Generation area 97

C

Calculate Score button 378 CanGrow property, reporting 478 CaptionPosition property, reporting 475 carbon-14 labeled compounds 552 Categorical Factors 68 Caution symbol, Analysis pane 119-120 Ceiling selection, mass defect 168, 224 Cell Selection Mode command 278 Cell Selection Mode shortcut menu command 586 Character spacing property, reporting 475 Charge parameter 146-147, 150, 508 charge state parameter 177, 180 structure editor 550 charged structures 540 Chart Type parameter 324 Check All Down-Regulated Points command 358 Check All Up-Regulated Points command 358 Check icon, structure checking tool 546 Check Point command 331 Check Structure message box 546 CheckAlignment property, reporting 478 CheckBox design item, reporting 472 Checked property filter, Result Filters view 267 Checked property, reporting 478 CheckSumEnabled property 478 ChemSpider Results table 416 Chromatogram Peaks table 384 Chromatograms view shortcut menu commands 302 working with 304 chromatographic peak apex, red line indicating 394 Clear All button, Result Filters view 274 Clear Compound Annotation command 279 Clear Frozen Content command 302 Close Sub-Report icon, reporting 467 Collapse All Column Headers command 278

collapsible left pane Chromatograms view 253 Mass Spectrum view 308 Options pane, Result Charts view 311 Collect Small Segments parameter, Pie Chart 325 Color legend, Scatter Chart view 330 color map, selecting 568 color palettes, standard 569 color schemes, selecting for a report template 445 ColumnCount property, reporting 480 ColumnDirection property, reporting 479 columns, application table changing the order 584 stacking 584 columns, report aligning 467, 469 transposing 444 ColumnSpacing property, reporting 480 communication tests 611 compatibility, software xxii Compound Annotation Editor dialog box 259 Compound Area Corrections view 367 compound class editor 538 Compound Class Matches table 429 Compound Class Scoring node 225 Compound Classes library, modifying 535 Compound Discoverer application window 7 Compound Editor dialog box for editing mass list compounds 524 shortcut menu 543 using 500 compounds labeled with carbon-14 552 Compounds table 401 compounds, library of expected 498 Configuration page 565 contacting us xxvi Continuous View icon, reporting 493 Control sample type 69, 112 Copy command 330 Copy Image command 301 Copy Points command 301 Copy with Headers command 278 copying Chromatograms view as a table of data points 301 Chromatograms view as an image 301 Copy to Clipboard dialog box 146 elemental composition of a compound 144 result tables 278 study factors 82, 104 cover page, report 465 CoverPage section, reporting 465

Create Analog Trace node connections 158 parameter descriptions 182 Create FISh Trace node, parameter descriptions 184 Create Mass Trace node, parameter descriptions 186 Create Pattern Trace node, parameter descriptions 187 Create Report icon 440 CrossSectionBox design item 473 CrossSectionLine design item 473 CSV files 5, 146, 278, 516 CTRL key, using 298 Culture property, reporting 480 Current Workflow Issues pane 120 Custom Filter Selection dialog box, features 598 Custom Isotope Ratio Pattern option 147 custom layout 258 Customize Report dialog box 442

D

data points, minimum number for peak detection 232 Data properties, reporting 480 Data Source parameter 326 Data Source, Scatter Chart 330 DataField property, reporting 481 date and time information, reporting 446 deactivating the software license 560 deactivation request file 561-562 dealkylation step, description 235 decimal precision for Is Equal To filter 267 default layout, result file 259 Define from Elemental Composition Formula option 146 Degradants folder 31 deleting adducts, library 503 compounds, library 499 fragments, compound class 538 ion definitions, library 507 mass lists 518 raw data files, effect of 99 study factors 82 transformations, library 511 workflow nodes in the Workflow Tree pane 141 workspace sections, report template 462 dendograms 369 Description box, working with 29 Descriptive Statistics view 347 Deselect All Visible Points command 331 Deselect Point command 331 design items, report template adding to a report template 469

Section Report pane 471 Design properties, reporting 483 DetailSection, reporting 465 Detect Compounds node 191 deuternopia, colormap for 569 diamond pattern tool 45 Differential Analysis view 352 Dimension Lines icon, reporting 468 direct match annotations 307 directional arrows layout guide tool 45 Metabolika pathways 531 Discriminate By check boxes 254 Distance Function parameter 375 Distribute Selection to All command 303 double bonds 203 Down Arrow, using 298 downloading documents xviii dragging input files to the Analysis pane 118 dragging views 45

E

E and L folder 31 Edit button, Analysis pane 133 Edit Compound Annotation command 279 Edit Sub-Report icon 458 editors Adduct Editor 503 biological replicate factor editor 81 categorical factor editor 79 Compound Annotation Editor 259 compound class editor 538 Compound Editor 500 Fragment Editor 539 Isotope Ratio Editor 142 numeric factor editor 80 Transformation Editor 512 Elemental Composition parameter 147 elements, changing 551 EMF files 291 Enable Column Fixing command 278 Enable Column Fixing shortcut menu command 585 Enable Row Grouping command 582 Enabled property, reporting 479 End property, reporting 483 Environmental folder 31 error bars 337 exclamation symbol 107 exclusion list 282 executing runs in parallel 565

Execution State column Analysis Results page 129 Job Queue page 128 Expand All Column Headers command 278 Expected Compounds library exporting to an XML file 499 features 499 importing from an XML file 499 initial 498 opening 499 Expected Compounds per File table 395 Expected Features table 397 Expected Finder node automatic peak width detection algorithm 174 connections 160 using 16-17, 19 Expected Formulas table 398 experimental variables 68 Export > As Mass List command 289–290 Export > Compounds to a New mzVault Library command 279 Export > Single Compound to an Existing mzVault Library command 279 Export As Xcalibur Inclusion/Exclusion List command 279 Export Image As command 301 Export Points As command 302 Export Spectra node connections 158 parameter descriptions 164 Export to Excel dialog box 281 Export to Existing Mass List dialog box 290 Export to New Mass List dialog box 289 Export to Text File command 278 exporting compounds to a new mzVault library 288 data to a CSV file 282 mass list for the TraceFinder application 286 reports 492 single compound to an existing mzVault library 287 Xcalibur inclusion/exclusion lists 282 Extract Sample Information From Sample Names dialog box 85

F

Factor Unit box, numeric factor 80 factory default result file layout 257 Factory Defaults command 314, 320, 323, 329 Features table 410 Field Chooser icon, using to show or hide columns 585 file path, input files 106 Files for Analysis area 118 Fill Gaps node connections 28, 159 parameter descriptions 193 Filter By check boxes 254 Filter By check boxes, Chromatograms view 253 Filter By Mass Defect node 167 Filter By Scan Event node 170 Filter Centroids node 173 filter set file saving 268 summary 276 Filter Summary page 276 Find dialog box, reporting 491 Find Expected Compounds node 230 Find icon, reporting 491 Find Missing Files command 109 firewalls 612 FISh Scoring node connections 160 FISh coverage score 36 using 232 FISh Scoring Queue view 368 FISh Trace Fragments table 386 Fit Page icon, reporting 493 Fit Width icon, reporting 493 Floor selection, mass defect 168, 224 flowchart, reporting 437 fold change 352 folder structure 70 Font property, reporting 476 Food Research folder 31 ForeColor property, reporting 476 Forensics folder 31 FormatString property, reporting 476 Forward connection, molecular networking 602 Fractional Mass selection, mass defect 168 Fragment Editor 539 Fragment Mode parameter 185 fragments, number of 537 FreeStyle data-visualization application 146 Freeze Content command, Chromatograms view 302 freeze pane 581

G

GdiCharSet property, reporting 476 GdiVerticalFont property, reporting 476 Generate Expected Compounds node connections 160 using 234 Generate Molecular Networks node 226

generating reports 439 Get from Clipboard button, Isotope Ratio Editor 145 Get from Composition Formula button, Isotope Ratio Editor 145 global settings 565 Go to Same Item in Main Table command 280 gray status rectangle 389-390 green background, table cell with 435 green circles, KEGG database 360 green status rectangle 389-390 Group By check boxes 254 Group By check boxes, Chromatograms view 253 Group By Sample check box 255 Group Compounds node 195 Group Expected Compounds node connections 160 parameter descriptions 239 grouping table rows 582 groups, sample 94 guide tool 45

Η

heat map 371 Height property, reporting 483 Help menu 11 Wizard 98 hidden result tables 257 table columns 585 Hide Grid icon, reporting 468 Hierarchical Cluster Analysis view 369 HighChem Fragmentation Options view 577 Highest Charge State parameter 177 high-mass compounds, predicted composition for 204 Histogram Chart displaying histograms 313 parameters 314 hydrogen-to-carbon ratio 203 Hyperlink property, reporting 481

I

Identification Only sample type 69 image files adding pictures to a report template 472 changing the logo image in a report template 453 saving data points to 291 selecting the logo image for a report template 445, 450 Image property, reporting 481 images of graphs 291

importing adduct libraries 503 compound library 5 ion definition libraries 507 library compounds 499 mzVault libraries 527 transformation libraries 511 Impurities folder 31 In ChemSpider column, Predicted Compositions table 207 inclusion list 282 inclusion/exclusion lists, exporting 282 indenting a related table in report templates 444 indenting columns in report templates 449 indicator icon, Result Filters window 269 Individual Traces parameter 184 Input File Characterization page 77 Input File Selection page 75 input files adding to a new study 75 dragging to the Analysis pane 118 modifying the location of 107 Input Files page of a study 105 Input Files table 387 installation instructions xxii Internet Explorer, set as default browser 308 Ion Definition Editor dialog box 509 Ion Definitions library adding ion definition 503 adding ion definitions 508 deleting entries 507 description 505 editing 509 exporting 507 features 507 importing ion definitions 503 Ion Definitions view 507 IP addresses, blocked 611 isolation mode, Molecular Networks viewer 606 Isotope check box 552 isotope pattern brackets, using to identify isotopic atoms 144 custom 147 intensity tolerance 191, 204 Pattern List Editor 149 reevaluating precursors 175 Isotope Ratio Editor, using 142 Isotope Ratios parameter 188 isotopologues 407 Isotopologues Distribution Chart 379

J

Job Queue page 124 jobs, processing in parallel 574

K

KeepTogether property, reporting 479 KEGG Pathways table 426 Kendrick Formula 168, 224 Known Structure Formats list 541

L

Label design item, reporting 471 Label Style parameter 324 labeled compounds 144, 188 Labeled Compounds per File table 414 Labeled Features table 412 Labeling Status column 407 Labels Font parameter 325 landscape paper orientation, report template 449 Last Changed parameter 104 layout files, result 257 Layout properties, reporting 483 Leaving Group parameter 514 result table column 400 Legend Size command 303 libraries adducts 502 compound classes 535 expected compounds 498 ion definitions 505 mass lists 515 transformations 510 license activation methods online 557 License Deactivation dialog box 560 License Manager page 553 license, software activation 558 deactivation 560 Life Sciences Mass Spectrometry Software Download and Licensing Portal 559 light bulb icon, New Study and Analysis Wizard 98 Line design item 472 Line Weight property, reporting 476 LineColor property, reporting 476 lines, adding to report templates 443 LineSpacing property, reporting 476 LineStyle property, reporting 476

Linkage Method parameter 375 Load button, Scatter Chart view 329 Load Filter Set dialog box 274 Location property, reporting 483 logo image changing 453 report templates 442 selecting 445 Lowest Charge State parameter 177

Μ

Manual Peak Integration command 303 Manual Peaks table 387 manual peaks, comparing to an XIC trace 22 Map to BioCyc Pathways node 217 Map to KEGG Pathways node 219 Map to Metabolika Pathways node 220 Mark Background Compounds node connections 28, 160 parameter descriptions 240 Mass Analyzer parameter 178 Mass Defect Plot view 339 Mass Defect Type parameter 168 Mass List Search Results table 418 mass lists exporting compounds to a new mass list 289-290 library 515 Mass Lists library 515, 520 Mass Lists view 518 mass spectrometers, HRAM 193 Max. Collision Energy parameter 171, 179 Max. Concurrent Entries, Xcalibur exclusion list 285 Max. Depth parameter 265 Max. Occurrence parameter 514 Max. Precursor Mass parameter 177 Max. Wavelength parameter 183 Maximum Shift parameter 166 MaxLength property, reporting 481 menu bar, application 7 Merge Features node 22, 241 Merged Features table 388 Metabolika Pathways table 427 Metabolika Pathways view 364 Metabolism folder 32 Metabolomics folder 32 Min. Collision Energy parameter 171, 179 Min. Precursor Mass parameter 177 Min. Wavelength parameter 183 Minimum Peak Count parameter 177 minimum peak intensity setting, recommended 193

Miscellaneous properties, reporting 484 missing files, resolving 107 mobile phase, background compounds in 240 MS Order parameter 170, 178, 186, 189 MS2 column 408 MultiLine property, reporting 479 Multipage view icon, reporting 493 multiple chromatograms, overlaying 298 mzCloud Results table 419 mzCloud search node 209 mzCloud Similarity button 378 mzCloud, submitting a single sample to 574 mzLogic Analysis view 377 mzLogic score 377 mzVault libraries adding compounds 287 creating new 288 mzVault Results table 421

Ν

n/a assignment, study factor 92 NarrowBarWidth property, reporting 476 nested experiments 68 New Analysis command, study 102 New Study and Analysis Wizard Extract Sample Information From Sample Names dialog box 85 Input File Characterization page 77 Input File Selection page 75 light bulb icon 98 Sample Groups and Ratios page 92 Study Name and Processing Workflow page 72 NewColumn property, reporting 481 NewPage property, reporting 482 nodes, workflow connections 158 parameters 163 Nominal Mass Rounding parameter 168, 224 normalization, batch 155 Normalize Areas node connections 28, 159 parameter descriptions 196 normalized areas, viewing 351 Nucleon Number list 552 Numerical Factors 68 NWRatio property, reporting 476

0

Open Analysis Template command, study 102 Open Common command, Workflows page 138 Open Response File dialog box 560 opening Atom Properties dialog box 551 existing studies 100 report templates 440 result files 248 structure files 541 Options panel, Scatter Chart view 330 OR logical conjunction, filtering with 271-272, 274 orange background, table cell with 435 orange status rectangle 389 Order, result table column 401 Orientation parameter, Custom Report dialog box 449 orientation, report template page 442 OutputFormat property, reporting 477 Overlay Cell Size box, Metabolika Pathways view 365 overlaying multiple chromatograms in one plot 298

Ρ

Padding property, reporting 483 page break, adding to a report 462 page numbers, adding to report templates 446 page orientation, selecting for a report 445 Page Thumbnails pane, reporting 495 PageBreak design item, reporting 472 PageFooter section, reporting 465 PageHeader bar, reporting 455 PageHeader section, reporting 465 Pan Mode icon, reporting 469, 494 paper size PaperKind parameter, Custom Report dialog box 442, 449 selecting for a report 445 Parallel Options view 565 parallel processing 574 Partial Least Squares - Discriminant Analysis view 358 Pattern List Editor 149 Pattern Scoring node connections 160 parameter descriptions 229 pausing processing runs 125 PDA detector, chromatograms 182 PDF file type, exporting reports 492 Phase parameter 514 result table column 400 Picture design item, reporting 472 PictureAlignment property, reporting 477 pie charts, parameter descriptions 323

pinning Options panel, Scatter Chart view 313 result table columns 278 study or result files on the Start Page 40 table rows 581 plot grid, Scatter Chart view 330 Polarity Mode parameter 172, 180 Polarity parameter 186, 189 polarity switching data 18 polarity, ionization 120, 184, 186, 189 portrait paper orientation, report template 449 Predict Compositions node 202 Predicted Compositions table 422 presentations, images for 291 Principal Component Analysis view 341 PrintAtBottom property, reporting 479 printing reports 492 Process Response File button 560 Processing Message Summary page 276 product ID for software license 555 Progress column, Job Queue 128 protanopia, colormap for 569 proxy server 614 Public Documents folder 65 publications, images for 291 p-values Differential Analysis node 26 volcano plot 352

0

QRCode property, reporting 485 Quality Control sample type 69 quickstart, Compound Discoverer 49 QuietZone property, reporting 477

R

Radius property, reporting 483 ratios, generating 97 RDBE calculation 203 red background, table cell with 435 red bar, heat map 370 red circle with exclamation symbol 107 red circles, KEGG database 360 red line, vertical 394 red status rectangle 389–390 Refresh button, Scatter Chart view 330 Refresh command 312 Region and Language settings xxi Regular Expression Builder area, Extract Sample Information from Sample Names dialog box 91 regular expression operators 88 Related Structures table 399 related tables basic targeted workflow 23 indenting in a report template 444 untargeted workflow 27 Relative Intensity command 303 Relative Mass Defect parameter selection 168 Remove All Plots command 302 Remove Analysis Result Files message box 107 Remove Files command, study 101 Remove Input File message box 107 Remove Plot command 302 removing chromatogram plots 297, 302 color schemes 447, 450 completed runs from the job queue 126 input files from the Input Files page 107 result or study files from the Start Page 40 RepeatToFill property, reporting 479 report designer page description 450 opening 453 shortcut menu 470 Report Preview dialog box, reporting 489 Report Print dialog box, reporting 489 report resolution page 491 report resolution page, reporting 489 ReportInfo design item, reporting 473 reporting Customize Report dialog box, using 442 exporting reports 492 previewing reports 489 printing reports 492 properties Appearance 474 Behavior 478 Data 480 Design 483 Lavout 483 Miscellaneous 484 report templates creating 442 data-and-time stamp, adding to reports 454 editing 450 page numbers 446 Reprocess command, Analysis Results page 129 reprocessing an analysis 131 request file, activation process 558 requirements

minimum number of data points for peak detection 232 software and hardware xix Reset button, Customize Report dialog box 450 resetting factory default layout, result file 257 Resolution parameter 147 restoring, factory default layout 257 Result Charts view 311 result files naming 133 opening 61, 248 Result Filters view 266 result tables BioCyc Pathways 217 BioCyc Results 425, 428 ChemSpider Results 416 Chromatogram Peaks 384 Compound Class Matches 429 Compounds 401 Expected Compounds per File 395 Expected Features 397 Expected Formulas 398 Features 410 FISh Trace Fragments 386 Input Files 387 KEGG Pathways 426 Labeled Compounds per File 414 Labeled Features 412 Manual Peaks 387 Mass List Search Results 418 Merged Features 388 Metabolika Pathways 427 Metabolika Results 428 mzCloud Results 419 mzVault Results 421 Predicted Compositions 422 Related Structures 399 Specialized Traces 390 Structure Proposals 391 Transformations 400 resuming processing runs 125 Retention Time Correction view 365 Reverse connection, molecular networking 602 reviewing analyses 129 result files 247 RichTextBox design item, reporting 472 RightToLeft property, reporting 479 ring equivalents 203 Rotation property, reporting 477 Round selection, mass defect 168, 224 Row Selection Mode, shortcut menu selection 144 rows, table copying 586 pinning 581 sorting 580 RT Offset parameter 182 Run command 122 Run command, Analysis pane 133 runs parallel processing 565 pausing 125 processing 125

S

S/N Threshold parameter 180 Sample Groups and Ratios page 92 sample groups and ratios, analysis summary 275 sample types 112 Samples page of a study 110 Save As icon, report designer page 466 Save Common command, Workflows page 138 Save Filter dialog box 274 saving all changes to open pages 8 analysis templates 133 auto-save option for studies 567 BioCyc credentials 576 current page 8 data points to a text file 5 data points to image files 5 filter sets 268 processing workflow templates 138, 161 report templates 466 result file layout 257 study factors 80, 82, 104 scaling factor, chromatographic peak area 199 Scan for Missing Features button, License Manager 563 Scan Polarity parameter 184 Scan Type parameter 172, 179 screen resolution, recommended xxi Scripting Node 244 Search ChemSpider node 206 Search Mass Lists node 207 Search mzCloud node 209 Search mzVault node 214 Search Results icon, reporting 495 sections, report designer page 464 Select All Visible Points command 331 Select Mode icon, reporting 468 Select Spectra node connections 158 parameter descriptions 174

using 174 Select Template Study File dialog box 74 Select/Deselect Item in All Groups icon 97 Selection Mode icon, reporting 492, 494 selection mode, choosing 549 Selection Tool, using 549 Set As Input File To shortcut menu command 112 Set Sample Type To command 85, 111 Shape design item, reporting 472 SHIFT key, using 298, 579 shifted match annotations 307 shortcut menu commands Chromatogram View 300 Compound Editor dialog box 543 Mass Spectrum View 310 report designer page 470 result tables 278 tabbed documents 44 Show All Tables check box, Result Filters view 267, 274 Show Details icon, Input Files page 106, 108 Show Detected Peaks command 303 Show Dots icon, reporting 468 Show Gridlines command 303 Show Legend command 303 Show Lines icon, reporting 468 Show Standard Errors command 337 Show Tooltips command 303 ShrinkToFit property, reporting 480 Single Page View icon, reporting 493 Size property, reporting 483 SizeMode, reporting 483 sizing bar, report designer page 462, 464 Sleeping status, Job Queue 127 Small Slice Collection parameter, Pie Chart 325 Small Slice Threshold (%) parameter 325 Snap Lines icon, reporting 468 Snap to Grid icon, reporting 468 Snapshot Mode icon, reporting 492, 494 software requirements xix sorting analyses on the Job Queue page 127 distribution maps, result table 580 result table rows 580 Sorting icon, study variable 97 Specialized Traces table 390 Specify FISh Scoring Settings dialog box 263 Spectrum Maximum parameter 183 spreadsheet exporting data to 278, 281 exporting the contents of a report 492

stable isotope labeling experiments 189, 379, 407 Stable Isotope Labeling folder 32 Standard Mass Defect parameter selection 168 Standard sample type 69, 112 Start Page hiding 41 icon 13 pin icon 40 recent file lists 40 shortcut menu 44 Start property, reporting 484 state, compound class file 538 status columns Expected Finder, Consolidated Peaks table 390 Unknown Detector, Consolidated Peaks table 389 status conditions, result filter 270 structure cleaning 548 creating a mirror image 547 moving 550 resizing 547 rotating 547 saving 550 structure drawing tools, using 542 Structure Editor module Atom Properties dialog box 550 bond multiplicity 550 bond properties 550 check structure 546 clean structure 548 mirror structure 547 move structure 550 resize structure 547 structure files, opening 541 Structure Proposals table 391 structures, charged 540 structures, checking the validity 546 studies Analysis Results page 129 Samples page 110 Studies Folder parameter 70 Study Name and Processing Workflow page 72 Study Summary page 276 study variables assigning 77 hierarchy 95 Style property, reporting 477 SubReport design item, editing 467 summaries, processing 275 Summed Trace parameter 184 SupplementOptions property, reporting 477 synchronizing the computer's clock 615

Т

table rows copying 586 grouping 582 pinning 581 sorting 580 tables filtering 588 freezing columns 585 sorting 580 Tag property, reporting 482 template tool, using 544 text file 5 Text property, reporting 482 Text with crossout line button, using 91 TextJustify property, reporting 477 Thermo MS Licensing, email from 555 Thermo Scientific website, user documents xviii Thumbnails Pane icon, reporting 495 time zone settings 615 Toggle Sidebar icon, reporting 493 tool, guide or position 45 toolbar, application 12 tooltips 336 Total Intensity Threshold parameter 177 Total Scan parameter 182 Trace Type parameter 182, 186 TraceFinder application, mass list for 286 Transformation Editor 512 transformations adding 512 editing 514 exporting 511 importing 511 Transformations library description 510 features 511 Transformations table 400 Transformations view 511 Transpose Data Fields shortcut menu command 470 transposing columns in a report template 469 transposing the columns in a report template 444 Trend Chart view 332 tritanopia, colormap for 569 troubleshooting access to the Online databases 611 analyses 57 Analysis pane Caution symbol 120 Current Workflow Issues pane 120 Region and Language settings xxi screen resolution xxi

U

Undo All Zoom/Pan command 300, 330 Undo Last Zoom/Pan command 300 units, chromatographic peak area 3 unrecognized scan event filters 180 Use Fragmentation Libraries check box 265 Use Isotope Pattern in Precursor Reevaluation parameter 175 Use Libraries parameter 185 Use Normalized Areas check box 351 UV-Vis detectors, chromatograms 182

V

validation issues missing Assign Compound Annotations node 123 missing Differential Analysis node 123 vector images 291 verbose messages, displaying in Job Queue 126 VerticalAlignment property, reporting 477 VerticalText property, reporting 477 View menu 9 viewer, molecular networking 604 views Bar Chart page, Result Charts view 319 BioCyc Pathways 362 Chromatograms 294 Compound Area Corrections 367 copying data to the Clipboard 291 Descriptive Statistics 347 FISh Scoring Queue 368 Hierarchical Cluster Analysis view 369 Histogram Chart page, Result Charts view 313 Isotopologues Distribution Chart 379 Job Queue page 124 KEGG Pathways 360 Mass Defect Plot view 339 Metabolika Pathways view 364 mzLogic Analysis view 377 Partial Least Squares - Discriminant Analysis 358 Pie Chart page, Result Charts view 323 Principal Component Analysis 341 Result Charts 311 Result Filters 266 Retention Time Correction 365 Scatter Chart page, Result Charts view 326 Trend Chart 332 Visible property, reporting 480

W

Wavelength Range parameter 183 whisker 333

width logo image container 450 object, report template 483 Overlay Cell Size box, Metabolika Pathways view 365 page width, report template 489 window position tool 45 workflow nodes installing 563 list of 163 numbers, hiding 567 parameters for Align Retention Times node 166 Analyze Labeled Compounds 189 Apply mzLogic node 223 Assign Compound Annotations node 200 Compound Class Scoring node 225 Create Analog Trace node 182 Create FISh Trace node 184 Create Mass Trace node 186 Create Pattern Trace node 187 Detect Compounds node 191 Export Spectra node 164 Export Xcalibur Inclusion/Exclusion List node 243 Fill Gaps node 193 Filter By Mass Defect node 167 Filter By Scan Event node 170 Filter Centroids node 173 Find Expected Compounds node 230 FISh Scoring node 232 Generate Expected Compounds node 234 Generate Molecular Networks node 226 Group Compounds node 195 Group Expected Compounds node 239 Map to BioCyc Pathways node 217 Map to KEGG Pathways node 219 Map to Metabolika Pathways node 220 Merge Features node 241 Normalize Areas node 196 Pattern Scoring node 229 Predict Compositions node 202 Scripting Node 244 Search ChemSpider node 206 Search Mass Lists node 207 Search mzCloud node 209 Search mzVault node 214 Select Spectra node 174 workspace sections, report designer page 464 Wrapmode property, reporting 477

X

X Axis Type parameter, Scatter Chart 329 X marks 302 Xcalibur Inclusion/Exclusion list, exporting 282

Y

Y Axis Type parameter, Scatter Chart 329 yellow background, table cell with 435 yellow-green background, table cell with 435

Ζ

Z Axis Type parameter, Scatter Chart 329 Zoom In icon, reporting 468 Zoom Out command 330 Zoom Out icon, reporting 468, 493 z-score transformation 374 Index: Z