

Thermo

TraceFinder

User Guide

Software Version 4.1
Optimized for Clinical Research

XCALI-97839 Revision A May 2016



Thermo
SCIENTIFIC

© 2016 Thermo Fisher Scientific Inc. All rights reserved.

TraceFinder, Aria, Q Exactive, FreeStyle, LCquan, ToxID, ExactFinder, Prelude, ISQ, and mzVault are trademarks, and Exactive, Thermo Scientific, TSQ Quantum, TSQ Endura, TSQ Quantiva, TurboFlow, and Xcalibur are registered trademarks of Thermo Fisher Scientific Inc. in the United States.

NIST is a registered trademark of the National Institute of Standards and Technology in the United States.

The following are registered trademarks in the United States and other countries: Windows, Active Directory, Excel, and Microsoft are registered trademarks of Microsoft Corporation. Waters and ACQUITY are registered trademarks of Waters Corporation.

ChemSpider is a trademark of ChemZoo, Inc.

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, May 2016

Software version: Microsoft Windows 7 Professional SP1; (Thermo) Foundation 3.1 SP3; Xcalibur 4.0; LC Devices 3.0; GC Devices 3.2

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

Contents

	Preface	vii
	Accessing Documentation	vii
	License Activation and Deactivation	ix
	Special Notices	x
	Contacting Us	xi
Chapter 1	Introduction	1
	About the TraceFinder Application	1
	TraceFinder Summary of Features	3
	TraceFinder Workflow	5
	Reporting Features	6
Chapter 2	Getting Started	9
	Installing the TraceFinder Application	9
	Installing the NIST and QED Libraries	14
	Launching the NIST Library Browser	17
	Launching a Qualitative Explorer	18
	Thermo Scientific FreeStyle	18
	Thermo Xcalibur Qual Browser	19
	Converting Legacy Data	20
	Converting Methods	22
	Converting Batches	24
	Converting Method Templates	26
	Converting Batch Templates	28
	Converting Compound Databases	30
	Choosing a Mode or Console	31
Chapter 3	Using the Acquisition Mode	35
	Working with Batches	36
	Opening and Navigating the Acquisition Mode	36
	Creating and Submitting Batches	38

	Real Time Status Pane	76
	Real Time Status – Acquisition Page	77
	Real Time Status – Instrument Page	78
	Real Time Status – Queues Page	79
	Instrument Controls Page	84
	Real-Time Trace Display	89
	Sample Types	91
Chapter 4	Using the Analysis Mode for Quantitation Batches	93
	Working in the Batch View for Quantitation Batches	94
	Samples Page for Quantitation Batches	94
	Auto Samples Page for Quantitation Batches	131
	Reference Sample Page for Quantitation Batches	133
	Threshold Samples Page for Quantitation Batches	134
	Working in Data Review for Quantitation Batches	135
	Sample View for Quantitation Batches	136
	Compound View for Quantitation Batches	151
	Compound Details for Quantitation Batches	164
	Comparative View for Quantitation Batches	192
	Qualitative View for Quantitation Batches	200
	Column Parameters for Compound Results and Sample Results	213
	Working in the Report View for Quantitation Batches	225
	Working in the Local Method View for Quantitation Batches	231
Chapter 5	Using the Analysis Mode for Target Screening Batches	233
	Working in the Batch View for Target Screening Batches	234
	Samples Page Features for Target Screening Batches	235
	Creating a New Target Screening Batch	249
	Editing a Target Screening Batch	256
	Submitting a Target Screening Batch	259
	Working in Data Review for Target Screening Batches	268
	Samples Pane	270
	Compounds Pane	271
	Chromatogram Pane	284
	Spectrum Pane	284
	Working in the Report View for Target Screening Batches	296
	Working in the Local Method View for Target Screening Batches	302
Chapter 6	Using the Analysis Mode for Unknown Screening Batches	305
	Working in the Batch View for Unknown Screening Batches	306
	Samples Page Features	306
	Creating a New Batch	318
	Editing a Batch	326
	Submitting a Batch	328
	Working in Data Review for Unknown Screening Batches	338

	Working in the Report View for Unknown Screening Batches	373
	Working in the Local Method View for Unknown Screening Batches	379
Chapter 7	Using the Report Designer.	381
	Editing a Template	382
	Toolbar Reference	394
	Home Toolbar	394
	Insert Toolbar – General Features	395
	Insert Toolbar – Quantitation Reports	396
	Insert Toolbar – Target Screening Reports	398
	Insert Toolbar – Unknown Screening Reports.	399
	Page Layout Toolbar	401
	Plot Display Properties	402
	Quantitation and Target Screening Plots.	402
	Unknown Screening Plots	418
	Quick Tips	430
Chapter 8	Using the Audit Viewer	431
	Audit Trail Log Files	431
	Audit Log File Functions.	432
	Audit Viewer	440
Appendix A	Using Quick Acquisition	443
Appendix B	Using Copy Down and Fill Down	445
Appendix C	Moving Data Review Panes	449

Preface

This guide describes how to use the Thermo TraceFinder™ 4.1 application with compatible Thermo Scientific™ GC/MS and LC/MS analytical instruments.

Contents

- [Accessing Documentation](#)
- [License Activation and Deactivation](#)
- [Special Notices](#)
- [Contacting Us](#)

❖ To suggest changes to the documentation or to the Help

Complete a brief survey about this document by clicking the button below.
Thank you in advance for your help.



Accessing Documentation

The TraceFinder application includes complete documentation. For system requirements, refer to the Release Notes on the software DVD.

❖ To view the TraceFinder manuals

From the Microsoft™ Windows™ taskbar, choose **Start > All Programs > Thermo TraceFinder > Manuals**.

—or—

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

❖ To view user documentation from the Thermo Fisher Scientific website

1. Go to thermofisher.com.
2. Click the **Services & Support** tab.
3. On the right, click **Manuals & Protocols**.
4. In the Refine Your Search box, search by the product name.
5. 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.

❖ To view TraceFinder Help

Open the TraceFinder application and choose **Help > TraceFinder Help**.

- To find a particular topic, use the Contents, Index, or Search panes.
- To create your own bookmarks, use the Favorites pane.

License Activation and Deactivation

Use the Thermo Scientific Product Licensing wizard to activate or deactivate the license for the TraceFinder application. To activate the license, you must have an activation code from Thermo Fisher Scientific. Before you transfer a license to another computer, deactivate the license.

When you first start the TraceFinder application, a dialog box displays the number of days remaining in your 120-day free evaluation license. If your evaluation license has expired, the License Activation wizard opens.

Note You can open the License Activation wizard at any time during your evaluation period by choosing **Help > About TraceFinder and Licensing** from the TraceFinder menu and then clicking **Activate**. If you already have a permanent license, a message tells you that your product is fully licensed.

Two types of licenses are available:

- 120-Day Evaluation Version (free of charge)
- Full Version Single License

The evaluation version is full-featured and automatically expires 120 days after activation. Any attempt to set back the system date automatically terminates this version. You can purchase and then activate the full version of the application at any time, during or after the free evaluation, without reinstalling the software.

Each activation key is valid only for a single license. Any additional installation generates a different license and requires a different activation key.

For software download and licensing questions, contact support at ThermoMSLicensing@thermo.com.

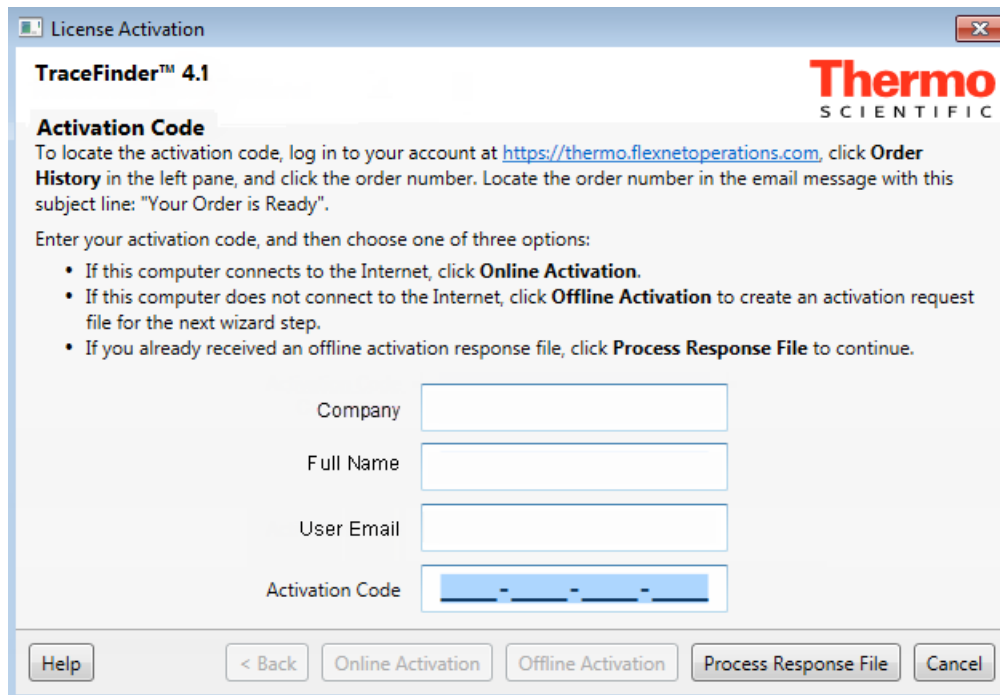
IMPORTANT The 120-day evaluation license includes the basic TraceFinder 4.1 features and the unknown screening features. When you purchase a permanent license, you have the option to purchase the unknown screening features. Your permanent license might not include the unknown screening features.

Use the License Activation wizard to activate or deactivate the license for the application. To activate the license, you must have an activation code from Thermo Fisher Scientific. You must deactivate the license before you transfer it to another computer.

❖ To start the license activation or deactivation process

1. Open the application.
2. Choose **Help > About TraceFinder and Licensing** to display the License Activation wizard.
3. Click **Activate (Deactivate)** to start the activation or deactivation process, as applicable.

The License Activation wizard opens.



4. Follow the instructions in the License Activation wizard.

For additional instructions, click **Help** in the wizard.

Special Notices

Make sure you follow the special notices presented in this guide. Special notices appear in boxes; those concerning safety or possible system damage also have corresponding caution symbols.

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

There are several ways to contact Thermo Fisher Scientific for the information you need. You can use your smartphone to scan a QR code, which opens your email application or browser.

Contact us	Customer Service and Sales	Technical Support
	(U.S.) 1 (800) 532-4752	(U.S.) 1 (800) 532-4752
	(U.S.) 1 (561) 688-8731	(U.S.) 1 (561) 688-8736
	us.customer-support.analyze@thermofisher.com 	us.techsupport.analyze@thermofisher.com 
 	<div><div>❖ To find global contact information or customize your request</div><div><div>1. Go to thermofisher.com.</div><div>2. Click Contact Us and then select the type of support you need.</div><div>3. At the prompt, type the product name.</div><div>4. Use the phone number or complete the online form.</div></div><div><div>❖ To find product support, knowledge bases, and resources</div><div>Go to thermofisher.com/us/en/home/technical-resources.</div></div><div><div>❖ To find product information</div><div>Go to thermofisher.com/us/en/home/brands/thermo-scientific.</div></div></div>	
<div><div>Note</div><div>To provide feedback for this document:</div><div><div>• Send an email message to Technical Publications (techpubs-lcms@thermofisher.com).</div><div>• Complete a survey at surveymonkey.com/s/PQM6P62.</div></div></div>		

Preface

Contacting Us

Introduction

This chapter describes general features of the TraceFinder application.

Contents

- [About the TraceFinder Application](#)
- [TraceFinder Summary of Features](#)
- [TraceFinder Workflow](#)
- [Reporting Features](#)

About the TraceFinder Application

The TraceFinder application targets the clinical research market. It supports a focused quantification workflow for specific nonbioanalytical laboratory use, instrument control, and method development functionality. TraceFinder is the primary application for the TSQ Quantum™ XLS triple quadrupole mass spectrometers.

The application can export mass data in the Acquisition List to XML format so that other systems' applications, such as for the TSQ 8000, TSQ Quantum™, ISQ, and Q Exactive™, can import the files into their databases.

The application can import the following file types:

- Sample lists in .csv or .xml format

See [Defining a Sample List](#) in [Chapter 3, "Using the Acquisition Mode."](#)

- Compounds from files that use the database (.xml or .cdb) format

Refer to Chapter 2, "Using Compound Databases in the Method Development Mode," in the *TraceFinder Lab Director User Guide*.

- Batches, methods, or templates from the TraceFinder 2.0, 2.1, 3.0, 3.1, 3.2, 3.3, or 4.0 applications.

See [Converting Legacy Data](#).

- Processing (.pmd) and instrument (.meth) method files from the Xcalibur™ data system

The *TraceFinder Lab Director User Guide* provides detailed information for creating the following methods.

Quantitative processing methods	Chapter 4, “Using the Method Development Mode for Quantitation Methods”
Target screening processing methods	Chapter 5, “Using the Method Development Mode for Target Screening Methods”
Unknown screening processing methods	Chapter 6, “Using the Method Development Mode for Unknown Screening Methods”
Instrument methods	Chapter 3, “Using Instrument Methods in the Method Development Mode”

The application checks the accuracy and precision of data against systems that have previously been certified against a standard processing program, such as the Statistical Analysis System (SAS).

Supported File Types

The application supports the following file types:

- Comma-separated values (.csv): A set of file formats used to store tabular data in which numbers and text are stored in plain textual form that can be read in a text editor. Lines in the text file represent rows of a table, and commas in a line separate fields in the tables row.
- Extensible Markup Language (.xml): A generic framework for storing any amount of text or any data whose structure can be represented as a tree. The only indispensable syntactical requirement is that the document has exactly one root element (also called the document element). This means that the text must be enclosed between a root start-tag and a corresponding end-tag.
- Instrument method (.meth): A proprietary file format for the Xcalibur data system suite with specific instructions that enable scientific instruments to perform data acquisition.
- Processing method (.pmd): A proprietary file format for the Xcalibur data system suite with specific instructions on processing data that was acquired through the instruments attached to the system.
- Raw data (.raw): The file type for acquired samples on the system.
- Compound database (.cdb): The file type for TraceFinder or ExactFinder™ compound database data.
- Library (.db): A library used for target screening.

TraceFinder Directory Structure

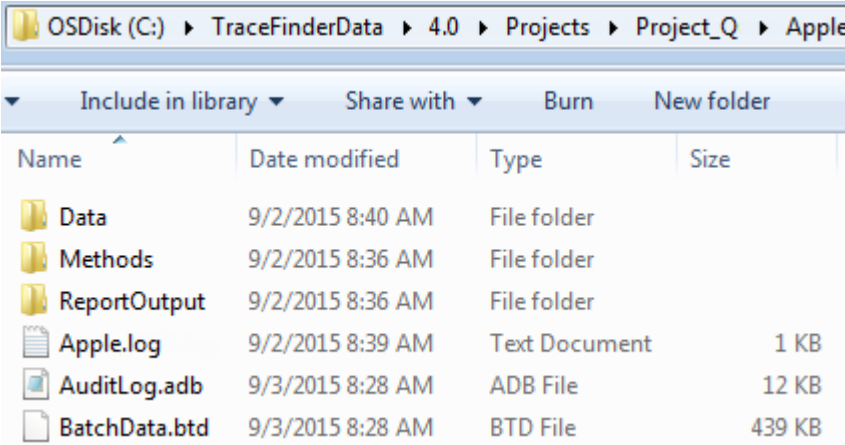
The application creates folders for batches, methods, and templates in the ...\\TraceFinderData directory. Within each batch folder, the application creates folders for data, methods, and reports.

IMPORTANT TraceFinder 4.1 uses the same data as TraceFinder 4.0. By default, the application stores the method, project, template, and compound database data for the 4.1 release in the TraceFinderData\\4.0 folder.

You can create batches in the TraceFinderData\\4.0\\Projects folder, or you can create subfolders within the Projects folder for your batches. You can create as many subfolders as you want for your batches, but you cannot create a batch within another batch folder.

IMPORTANT You cannot rename or move the folders created by the TraceFinder application.

Figure 1. Example batch directory structure



Name	Date modified	Type	Size
Data	9/2/2015 8:40 AM	File folder	
Methods	9/2/2015 8:36 AM	File folder	
ReportOutput	9/2/2015 8:36 AM	File folder	
Apple.log	9/2/2015 8:39 AM	Text Document	1 KB
AuditLog.adb	9/3/2015 8:28 AM	ADB File	12 KB
BatchData.btd	9/3/2015 8:28 AM	BTD File	439 KB

TraceFinder Summary of Features

The TraceFinder system provides a workflow-oriented approach to high-throughput quantitation, using tools that automate and speed up the processes of method creation, loading samples, automatically generating data, manually reviewing and editing results, and finalizing the data review and reporting process.

The TraceFinder software package includes data acquisition, processing, reviewing, and reporting capabilities designed to assist analysts in clinical research applications. The application has a fully automated acquisition mode and a manual data analysis mode. You can use the data acquisition system to create and submit batches and monitor real-time review of results.

The application uses a comprehensive processing method to provide improved handling of ion ratio calculations, reviewing, and reporting. In addition, it can compare the mass spectra and integrate the processes of data review and reporting.

Key features include the following:

- Role-based authorization for Security, LabDirector, ITAdmin, Supervisor, Technician, and QAQC (quality assurance) roles
- Administrator Console for user security, role-based permissions, and data repositories
- Configuration console for report configuration, detection and acquisition defaults, adduct definitions, screening library selection, and customized columns and flags
- Method Development mode for editing instrument methods, setting processing and error flag parameters, and setting reporting options
- Acquisition mode that guides you in creating batches and running samples
- Analysis mode with batch views, data review, local method views, and reporting views
- Database-capable method development
- Quantification, target screening, and unknown screening workflows
- Spreadsheet-based report designer

Features of the common workflow core include the following:

- Acquisition and processing
- Peak detection
- Quantification to include calibration
- Error analysis and flag setting
- Reporting
- Data persistence
- Raw data file handling

TraceFinder Workflow

The application is structured with a typical laboratory workflow in mind. You create a batch, and the system injects samples into the instrument, runs the samples, analyzes the data, and generates a report. You can set up a master method for specific compound groups or assays that you expect to run in your laboratory. When you are ready to run a particular type of sample, select the appropriate method and begin.

When using the application, follow these basic steps:

1. Create and save a master method in the Method Development mode.

A master method combines the instrument method and processing method that define the following:

- How the raw data is acquired and processed
- How the error checking information evaluates the results
- How the results appear in reports



2. Create and submit a batch using the Acquisition wizard.

A batch lists samples for processing and reporting using a specified method. Each row of a batch represents a unique sample.



3. Monitor the status of the batch in the Real Time Status view.

The real-time display is visible from all the TraceFinder modes. You can begin another batch while you watch the real-time display of the currently acquiring batch.

Note At any time, you can quickly view the system status by looking in the upper right corner of the TraceFinder window. This area displays a green, yellow, or red status light and a description of any activity in the queues, as in this example:

Apple_0_81_01  Acquiring 

The status light (in any color) might include an exclamation mark.

- An exclamation mark on a green status light, , might indicate that the instrument method contains the wrong source (for example, APCI instead of HESI).
- An exclamation mark on a red status light, , might indicate that a vial is missing and the batch was forced to stop.

Click the status light to display details of the error.

4. Evaluate the data in the Analysis mode.

The Analysis mode includes views where you can review batches, batch data, reports, and local methods.

5. View and print reports in the Report View of the Analysis mode.

Use the Report View to view or print the reports for the current batch.

Reporting Features

The report engine can generate several different types of reports designed to meet the needs of the laboratory, the laboratory's customers, and key regulatory agencies that might review the results. The following types of reports meet the requirements of various methods and global regulatory agencies, helping to track the performance of LC and GC systems and methods.

Standard Report Types

- Ad Hoc Tune Report
- Batch Report
- Blank Report
- Breakdown Report
- Calibration Report
- Check Standard Report
- Chromatogram Report
- Compound Calibration Report
- Compound Calibration Report - Alternate
- Confirmation Report
- Confirmation Report 2
- High Density Calibration Report
- High Density Internal Standard Report Long
- High Density Sample Report 1 Long
- Intelligent Sequencing
- Internal Standard Summary Report
- Ion Ratio Failure Report
- Manual Integration Report
- Method Report
- Method Validation Report
- MSMSD Report
- Quantitation Report
- Quantitation Report - 2
- Sample Report

- Solvent Blank Report
- Standard Addition Report
- Surrogate Recovery Report
- Target Screening Height Density Sample Report
- Target Screening Height Density Sample Report 2
- Target Screening Summary Report
- TIC Report
- TIC Summary Report
- Tune Report
- Unknown Screening High Density Report
- Unknown Screening Long Report
- Unknown Screening Summary Report

Use the features in the Report Designer to create custom report types. See [Chapter 7, “Using the Report Designer.”](#)

Getting Started

This chapter includes the procedures for getting started with the TraceFinder application.

Contents

- [Installing the TraceFinder Application](#)
- [Installing the NIST and QED Libraries](#)
- [Launching the NIST Library Browser](#)
- [Launching a Qualitative Explorer](#)
- [Converting Legacy Data](#)
- [Choosing a Mode or Console](#)

Installing the TraceFinder Application

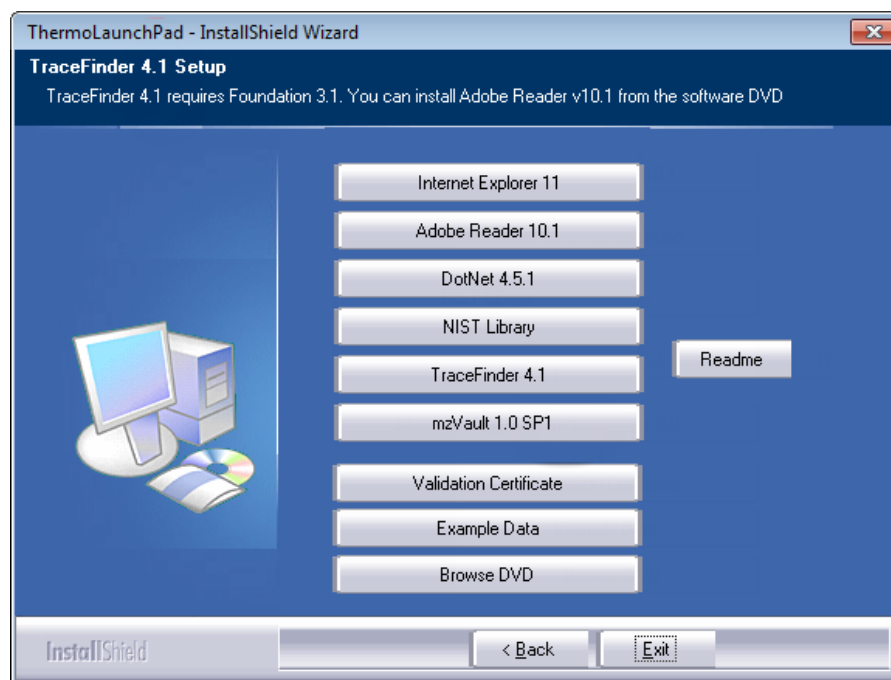
To initially install the TraceFinder 4.1 application, follow the instructions in the *TraceFinder Installation Guide*. Later, you might need to reinstall the application or other features using the InstallShield Wizard.

Follow these instructions to reinstall, start, and log in to the application.

❖ To reinstall the TraceFinder application

1. From the Thermo Foundation Instrument Configuration window, remove all instruments.
2. From the Windows Control Panel, uninstall the application and then uninstall all Thermo instrument drivers.
3. Insert the TraceFinder DVD, and install both the TraceFinder 4.1 application and the NIST™ library as follows:
 - a. Open the TraceFinder launcher and click **Next**.

The InstallShield Wizard opens.



- b. Click **TraceFinder 4.1**, and follow the instructions in the InstallShield Wizard.

The installer verifies that you have the appropriate versions of the Thermo Foundation™ platform and Xcalibur data system and updates them if necessary.

- c. At the prompt, click **Yes** to completely remove any previously installed applications.
- d. Open the TraceFinder launcher again and click **Next**.
- e. Click **TraceFinder 4.1**, and follow the instructions in the InstallShield Wizard.

IMPORTANT For the application to properly install, you might be prompted to uninstall Foundation platform. Do the following:

1. Click **Yes**, and then when prompted to restart your computer, click **OK**.

The wizard continues the TraceFinder installation.

2. When prompted to install Thermo Foundation, click **Yes**, and then when prompted to restart your computer, click **OK**.

The wizard continues the installation.

- f. When prompted, choose to install either the **GC** or **LC** version of the software.
- g. When the installation is complete, open the TraceFinder launcher again and click **Next**.
- h. When prompted to restart your computer, select the **Yes** option and click **Finish**.

3. (Optional) After your system restarts, open the TraceFinder launcher again, click **Next**, and install any of the following:
 - a. Click **NIST Library** and follow the instructions to install it.
 - b. Click **Example Data**.

The application installs example compound databases, instrument methods, and batch data.
 - c. Click **mzVault 1.0 SP1** and follow the instructions to install the Thermo mzVault application.
4. Install the appropriate device drivers, and configure the instruments in the Thermo Foundation Instrument Configuration window.

You can now start the TraceFinder application.

❖ **To start the TraceFinder application**

1. Configure your instruments.

You must close the TraceFinder application before you can configure your instruments.

2. Double-click the **TraceFinder** icon on your desktop, or choose **Start > All Programs > Thermo TraceFinder 4.1 > TraceFinder 4.1 Clinical Research**.

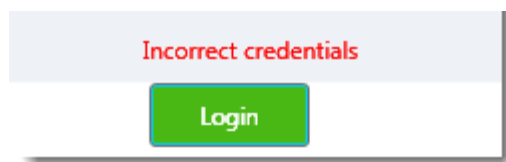
By default, user security is not activated and the application does not require a password. To activate user security, refer to the *TraceFinder Administrator Console User Guide*.

❖ **To log in to the TraceFinder application (when user security is activated)**

Note Before you can log in to the application when user security is activated, a system administrator must set up a user account for you.

1. Enter your user name in the [TraceFinder login window](#).
2. Enter your password.

If your user name or password does not match, the system reports this error:



Correct the user name or password, or contact your system administrator.

3. Click **Login**.

The [TraceFinder main window](#) opens.

Figure 2. TraceFinder main window, showing the Analysis mode for a quantitation batch

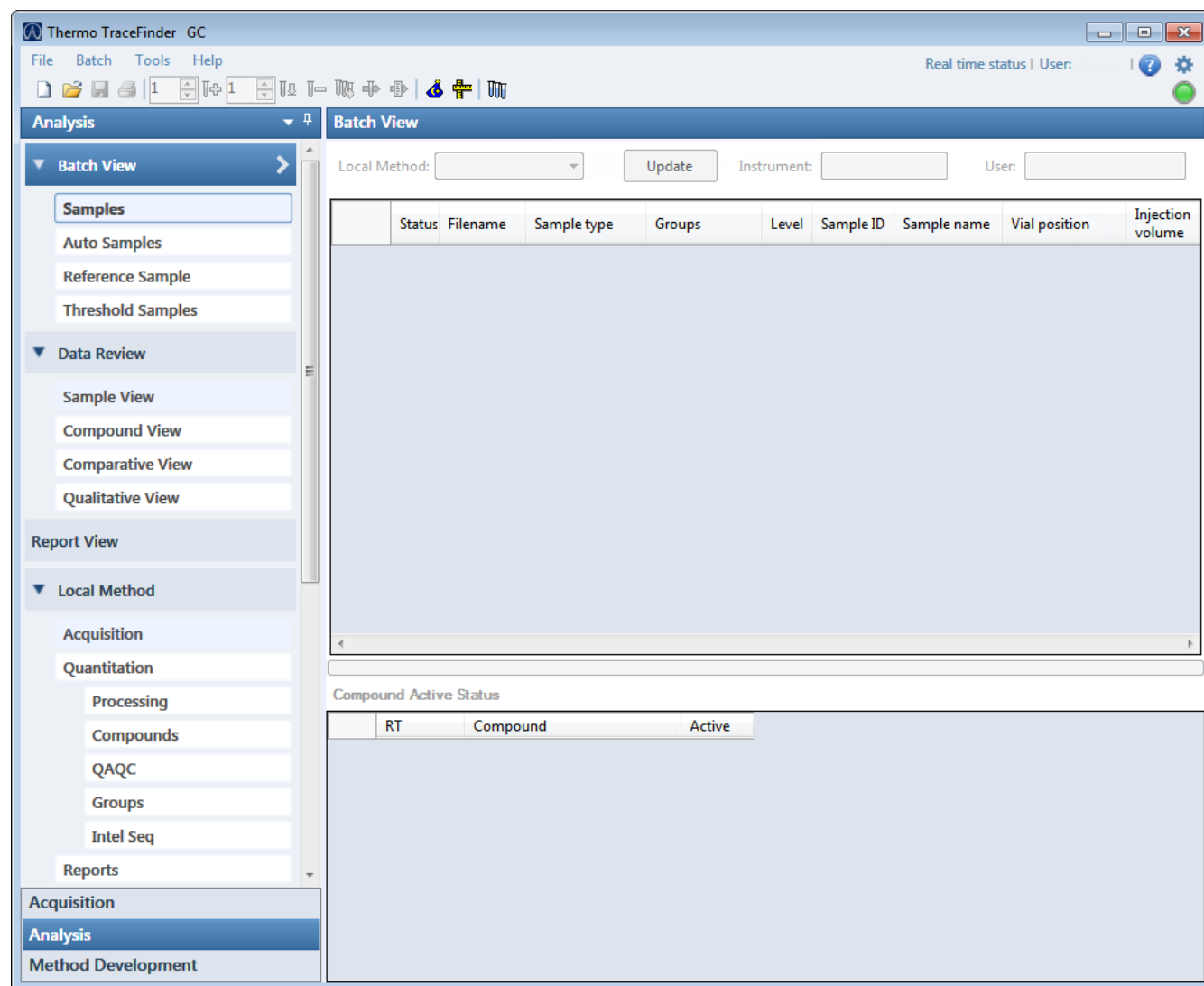


Table 1. TraceFinder main window features (Sheet 1 of 2)

Parameter	Description
Modes	
Acquisition	See Chapter 3, “Using the Acquisition Mode.”
Analysis	See Chapter 4, “Using the Analysis Mode for Quantitation Batches.”
	See Chapter 5, “Using the Analysis Mode for Target Screening Batches.”
	See Chapter 6, “Using the Analysis Mode for Unknown Screening Batches.”
Method Development	Refer to the <i>TraceFinder Lab Director User Guide</i> .

Table 1. TraceFinder main window features (Sheet 2 of 2)

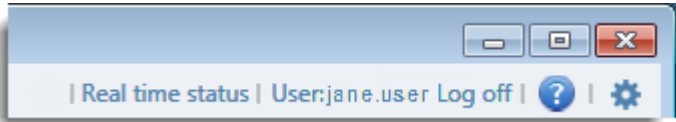

Parameter	Description
	A TraceFinder window with user security for a user in the default LabDirector role has these functions.
Additional Features	
Real Time Status	Opens the Real Time Status pane for the current acquisition. The acquisition progress is displayed within the current mode window.
User	Displays the name of the current user.
Log Off	Logs off the current user and displays the login screen. This function is available only when user security is activated.
Help ?	Opens the TraceFinder Help.
Application Configuration 	Opens the Configuration console where you can configure several options for using the application. Refer to Chapter 1, “Using the Configuration Console,” in the <i>TraceFinder Lab Director User Guide</i> .

Figure 3. TraceFinder login window

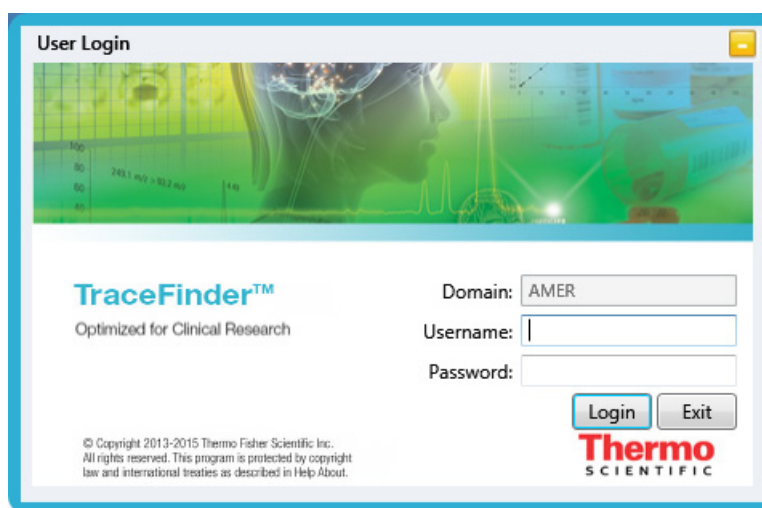


Table 2. Login window parameters

Parameter	Description
Domain	The authentication method.
Username	The user's assigned user name.
Password	The assigned password for the user name.
Login	Verifies the user name and password, and opens the application.
Exit	Closes the TraceFinder login window.

Installing the NIST and QED Libraries

When you are using triple quadrupole instruments, such as the TSQ Quantum XLS, follow these instructions to install the NIST and QED libraries.

❖ To install the NIST library


1. Open the TraceFinder launcher, and click **Next**.
2. Click **NIST Library**.

The NIST 08 MS Search and AMDIS Setup wizard opens.

3. Follow the instructions in the setup wizard.
4. When the wizard prompts you to select a destination folder, select **C:\Program Files\NISTMS**.
5. Continue to follow the instructions in the wizard until the setup is complete.

Note The application installs the NIST 2008 library; however, it also supports the NIST 2011 and NIST 2014 libraries.

❖ To install the QED library

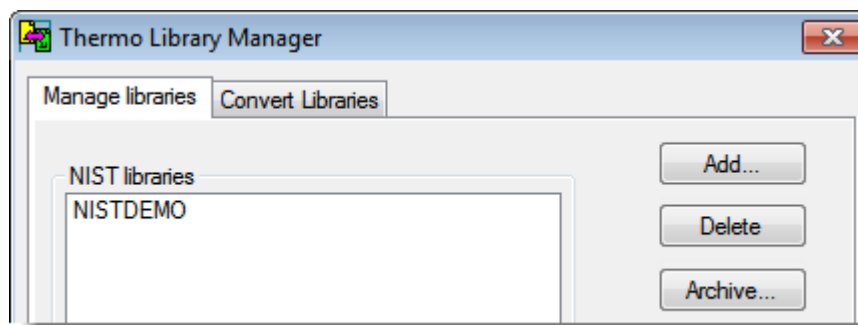
1. On your desktop, double-click the **Xcalibur** icon, .

The Thermo Xcalibur Roadmap opens.



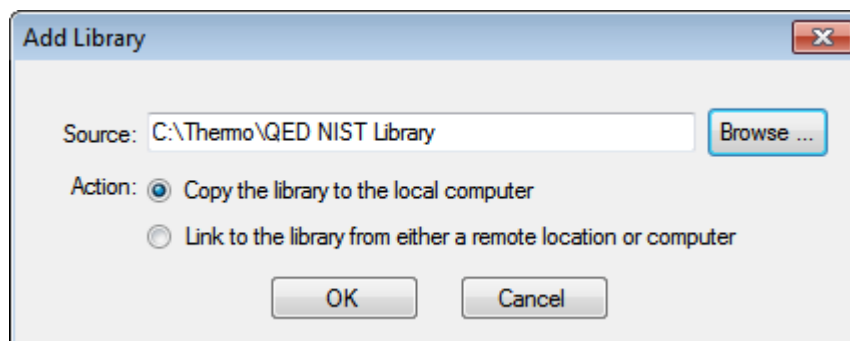
2. Choose **Tools > Library Manager** from the main menu.

The Thermo Library Manager dialog box opens, showing the NIST Libraries list.



3. Click **Add**.

The Add Library dialog box opens.

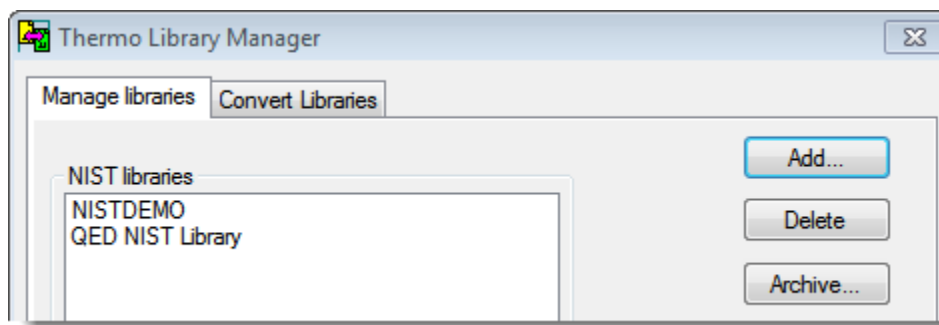


4. Click **Browse**, and locate your QED library in the C:\Thermo folder.
5. Click **OK**.

The Xcalibur data system reports that it has added the library to the NIST application.

6. Click **Dismiss** to close the message box.

The Xcalibur data system adds the QED library to the NIST Libraries list in the Library Manager dialog box.

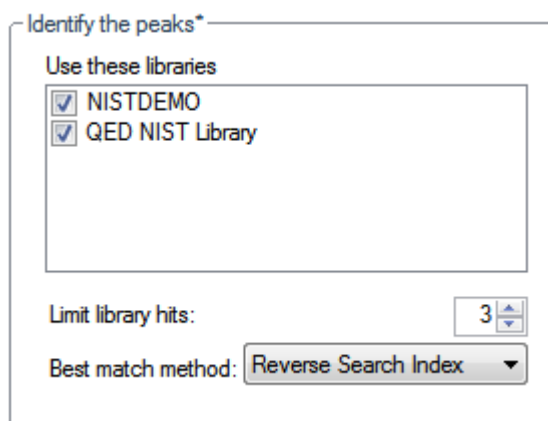


2 Getting Started

Installing the NIST and QED Libraries

7. Click **Exit** in the Thermo Library Manager dialog box.
8. To confirm the library installation, do the following:
 - a. Start the application.
 - b. Click **Method Development** in the navigation pane.
 - c. Click **Method View** in the Method Development navigation pane.
 - d. Choose **File > New > Method Template** from the main menu.

The Method Template Editor displays the QED NIST Library in the Use These Libraries list.

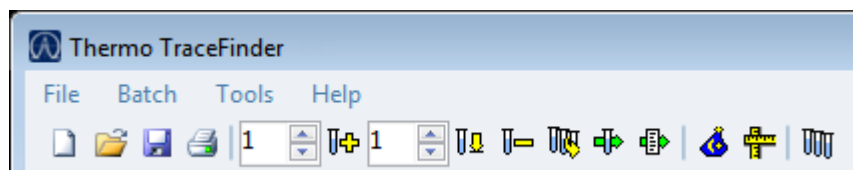


Launching the NIST Library Browser

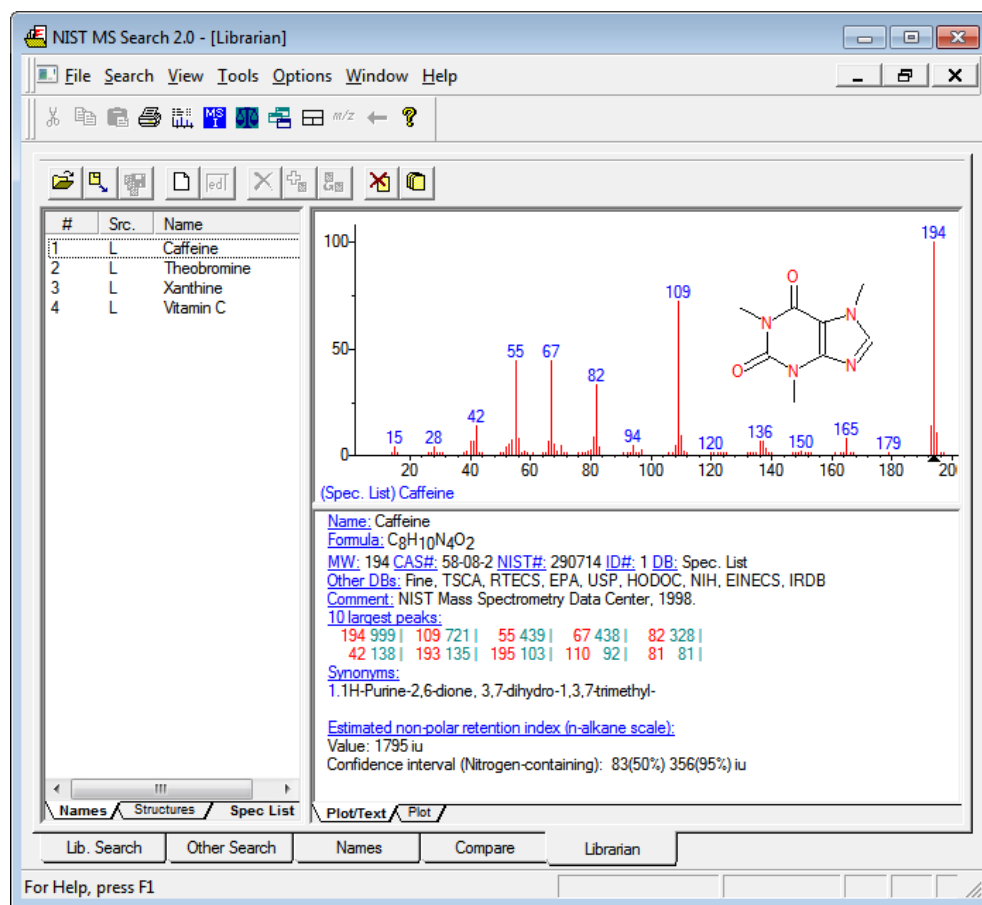
Use the NIST MS Search tool to search the NIST library.

❖ To open the NIST library browser

Choose **Tools > Launch Library Browser** from the TraceFinder main menu.



The NIST MS Search window opens.



For detailed instructions about using the library browser, refer to the Help in the NIST MS Search window.

Launching a Qualitative Explorer

Use a qualitative explorer application to display chromatograms and spectra, detect chromatogram peaks, search libraries, simulate spectra, subtract background spectra, apply filters, add text and graphics, create and save layouts, and view instrument parameters as they changed during the acquisition.

Your TraceFinder application is configured to use one of the following applications:

- Thermo Scientific FreeStyle
- Thermo Xcalibur Qual Browser

Thermo Scientific FreeStyle

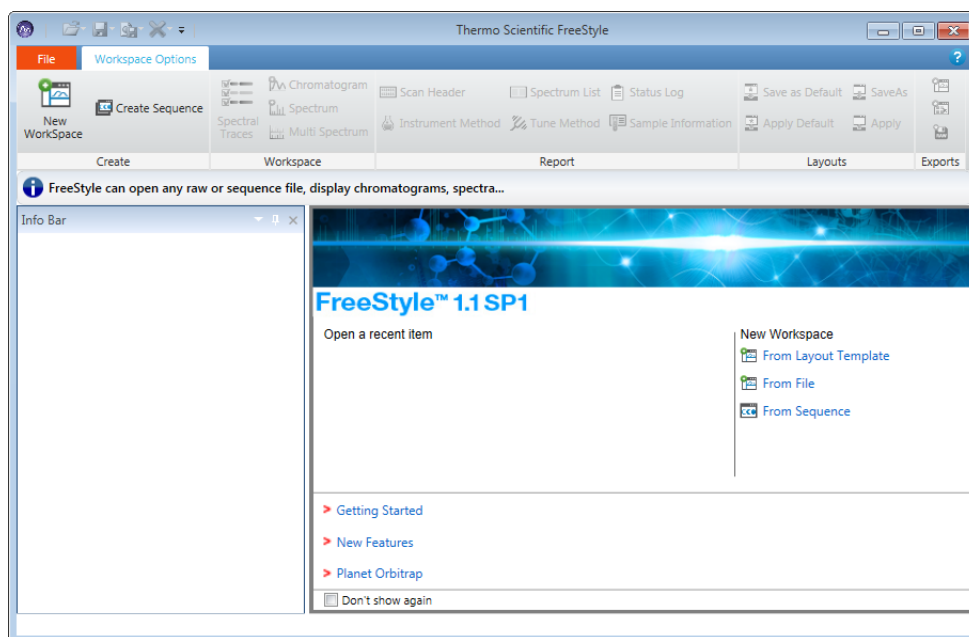
IMPORTANT The Thermo Scientific FreeStyle™ application is available only when you configure it as your default qualitative explorer in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.

❖ To open the FreeStyle window

Choose **Tools > Launch Qual Explorer** from the TraceFinder main menu.

The FreeStyle application opens.

Figure 4. FreeStyle main window



For detailed instructions about using the FreeStyle application, click the **Help** icon, , in the FreeStyle window.

Thermo Xcalibur Qual Browser

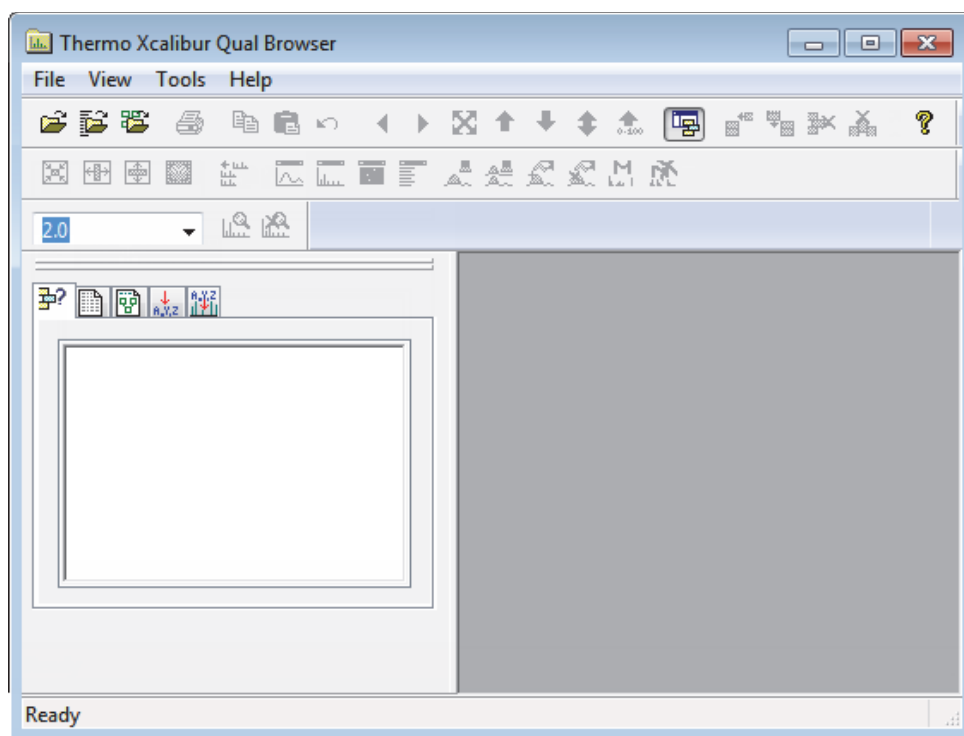
IMPORTANT The Qual Browser application is available only when you configure it as your default qualitative explorer in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.

❖ **To open the Qual Browser window**

Choose **Tools > Launch Qual Explorer** from the TraceFinder main menu.

The Thermo Xcalibur Qual Browser application opens.

Figure 5. Qual Browser main window



For detailed instructions about using the Qual Browser application, refer to the Help in the Qual Browser window.

Converting Legacy Data

Use the TraceFinder Legacy Data Converter to convert methods, batches, method templates, batch templates, or compound databases (CDBs) from the source versions to compatible TraceFinder 4.0/4.1 target configurations.

IMPORTANT TraceFinder 4.1 uses the same data configuration as TraceFinder 4.0. You do not need to convert your data from 4.0 to 4.1. Data for the 4.1 release is stored in the TraceFinderData/4.0 folder.

- You can convert legacy methods, batches, or method templates from TraceFinder versions 2.0, 2.1, 3.0, 3.1, 3.2, or 3.3.
- You can convert legacy batch templates from TraceFinder versions 3.0, 3.1, 3.2, or 3.3.
- You can convert legacy compound databases from TraceFinder versions 3.0, 3.1, 3.2, or 3.3.
- You can convert data from TraceFinder version 4.0/4.1 for general quantitation to another installed configuration of TraceFinder 4.0/4.1.

Version Compatibility

This table shows which source versions of methods, batches, method templates, batch templates, or compound databases are compatible with TraceFinder 4.0/4.1 target configurations.

Table 3. Version compatibility (Sheet 1 of 2)

Source	TraceFinder 4.0/4.1 target			
	General	EFS ^a	Clinical Research	Forensic Toxicology
TraceFinder 4.0/4.1 General		✓	✓	✓
TraceFinder 3.3 General		✓	✓	✓
TraceFinder 3.3 EFS		✓		
TraceFinder 3.3 Clinical Research			✓	✓
TraceFinder 3.3 Forensic Toxicology			✓	✓
TraceFinder 3.2 General		✓	✓	✓
TraceFinder 3.2 EFS		✓		
TraceFinder 3.2 Clinical Research			✓	✓
TraceFinder 3.2 Forensic Toxicology			✓	✓
TraceFinder 3.1 General		✓	✓	✓
TraceFinder 3.1 EFS		✓		
TraceFinder 3.1 Clinical Research			✓	✓
TraceFinder 3.1 Forensic Toxicology			✓	✓
TraceFinder 3.0 General	✓	✓	✓	✓

Table 3. Version compatibility (Sheet 2 of 2)

Source	TraceFinder 4.0/4.1 target			
	General	EFS ^a	Clinical Research	Forensic Toxicology
TraceFinder 3.0 EFS		✓		
TraceFinder 3.0 Clinical Research			✓	✓
TraceFinder 3.0 Forensic Toxicology			✓	✓
TraceFinder 2.1 General	✓	✓	✓	✓
TraceFinder 2.1 EFS		✓		
TraceFinder 2.1 Clinical Research			✓	✓
TraceFinder 2.1 Forensic Toxicology			✓	✓
TraceFinder 2.0 General	✓	✓	✓	✓
TraceFinder 2.0 EFS		✓		
TraceFinder 2.0 Clinical Research			✓	✓

^a Environmental and Food Safety

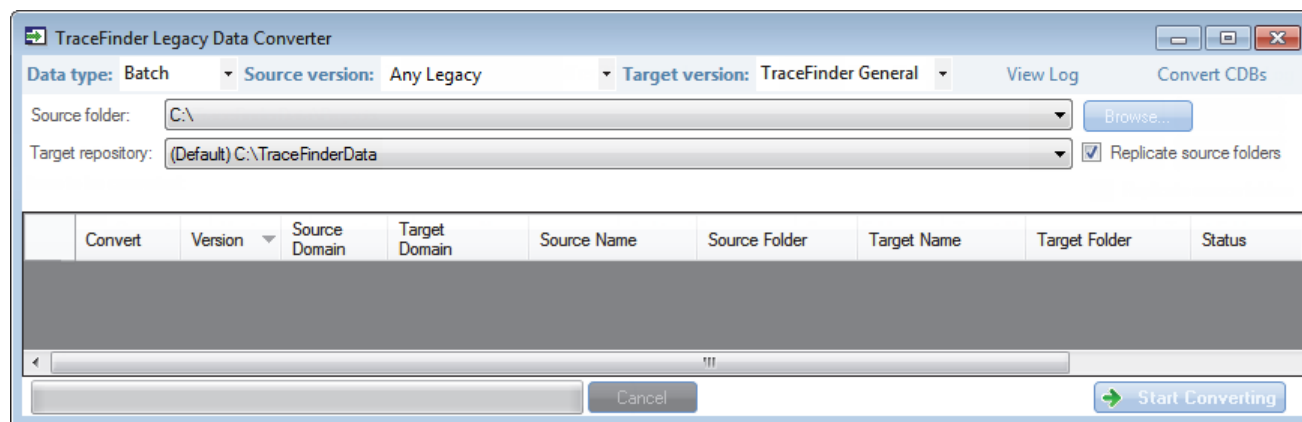
See the following topics:

- [Converting Methods](#)
- [Converting Batches](#)
- [Converting Method Templates](#)
- [Converting Batch Templates](#)
- [Converting Compound Databases](#)

❖ **To open the TraceFinder Legacy Data Converter**

Choose **Tools > Launch Legacy Data Converter** from the TraceFinder main menu.

The TraceFinder Legacy Data Converter window opens.



Note When you open the application, the system checks for any legacy data and prompts you to open the Legacy Data Converter.

Converting Methods

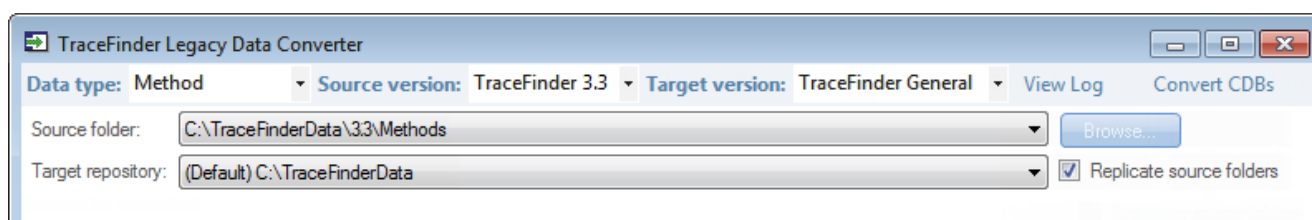
Use the data converter to convert legacy methods to TraceFinder 4.1 methods.

❖ To convert a method

1. In the Data Type list, select **Method**.

The TraceFinder Legacy Data Converter displays the interface for converting methods.

The following example shows that you can convert methods from the TraceFinder 3.3 General configuration to the current General configuration. For a complete list of version compatibilities, see [Version Compatibility](#).



2. In the Source Version list, select the version of the method that you will convert.

Note When you select Any Legacy, the Legacy Data Converter examines all possible methods in the source folder, regardless of version.

The conversion table displays the methods in the Methods folder for the selected source version. The application verifies that the method file is in the .mmx file format.

3. To convert a method that is not in the default list, do the following:
 - a. Click **Browse** and locate a different source method folder.

You can select a specific method folder or a folder that contains multiple methods.

- b. Click **OK** in the Browse for Folder dialog box.

The application displays the selected method folder in the conversion table.

When you select a folder that contains multiple method folders, the application displays all the methods.

4. In the Target Version list, select the version that you are converting to.

The list displays only TraceFinder configurations with compatible data. See [Version Compatibility](#).

5. (Optional) In the Target Name column, change the default new name for each method that you want converted.

When you populate the conversion table, the application checks each method to see if a method with this name exists in the target repository.

- If the method name already exists in the target repository, the default new name appends “_1” to the original name.
- If the method name does not exist in the target repository, the application keeps the original method name.

IMPORTANT The conversion cannot overwrite an existing file name. If the new name is identical to an existing method file, the conversion will not work. When you manually enter a new name, you must verify that the name does not already exist.

6. Select the **Convert** check box for each method that you will convert, and click



The application confirms that all methods to be converted use the .mmx file format.

When the conversion process begins, the application displays a status bar and a Cancel button. You can cancel pending conversions, but not the method that is currently converting.

When the Status column reports that a method is successfully converted, the application writes the converted file to the specified target repository.

Note If a method conversion is unsuccessful, the Status column displays “Conversion failed.” The log file contains details about the failed conversion.

7. To view a log of the conversion, click **View Log**.

The application opens a cumulative log file for the session in a Microsoft Notepad text editor window.

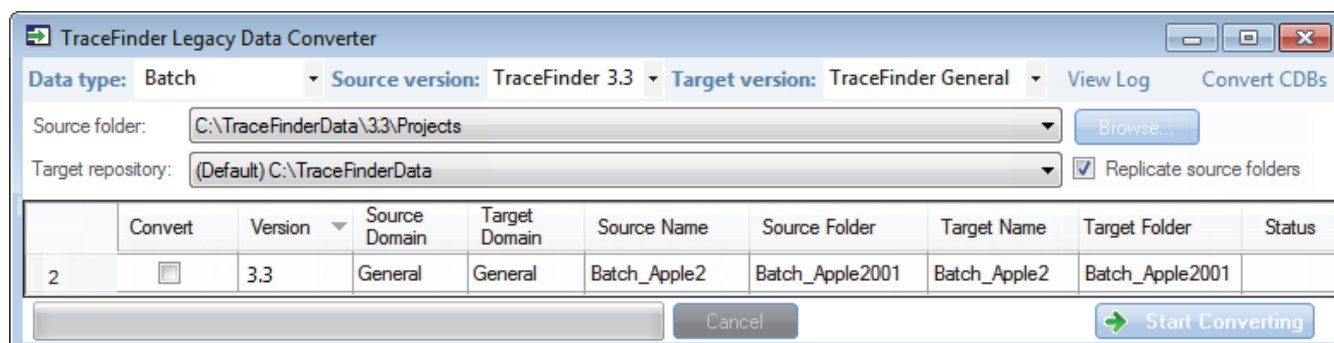
Converting Batches

Use the data converter to convert legacy batches to TraceFinder 4.1 batches.

❖ To convert a batch

1. In the Data Type list, select **Batch**.

The following example shows that you can convert batches from the TraceFinder 3.3 General configuration to the current General configuration. For a complete list of version compatibilities, see [Version Compatibility](#).



2. In the Source Version list, select the version of the batch that you will convert.

Note When you select Any Legacy, the Legacy Data Converter examines all possible batches in the source folder, regardless of version.

The conversion table displays all batches in the Projects folder for the selected source version.

3. In the Target Version list, select the version that you are converting to.

The list displays only TraceFinder configurations with compatible data. See [Version Compatibility](#).

4. In the Target Default Project and Subproject boxes, type the name of a project and subproject, or select the **Replicate Original Project/Subproject** check box.
5. (Optional) In the New Name column, change the default new name for each batch that you want converted.

When you populate the conversion table, the application checks each batch to see if a batch with this name exists in the target repository.

- If the batch name already exists in the target repository, the default new name appends “_1” to the original name.
- If the batch name does not exist in the target repository, the application keeps the original batch name.

IMPORTANT The conversion cannot overwrite an existing file name. If the new name is identical to an existing batch folder, the conversion will not work. When you manually enter a new name, you must verify that the name does not already exist.

6. Select the **Convert** check box for each batch that you will convert, and click



The application confirms that all batches to be converted use the .btx file format.

When the conversion process begins, the application displays a status bar and a Cancel button. You can cancel pending conversions, but not the batch that is currently converting.

When the Status column reports that a batch is successfully converted, the application writes the converted batch to the ...\\TraceFinderData\\4.0\\Projects folder and uses either the original project and subproject names or the new names that you entered.

Note If a batch conversion is unsuccessful, the Status column displays “Conversion failed.” The log file contains details about the failed conversion.

7. To view a log of the conversion, click **View Log**.

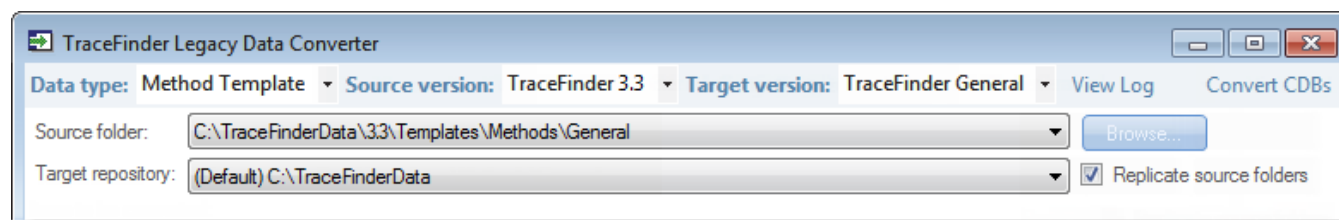
Converting Method Templates

Use the data converter to convert legacy method templates to TraceFinder 4.1 method templates.

❖ To convert a method template

1. In the Data Type list, select **Method Template**.

The following example shows that you can convert method templates from the TraceFinder 3.3 General configuration to the current General configuration. For a complete list of version compatibilities, see [Version Compatibility](#).



2. In the Source Version list, select the version of the method template that you will convert.

Note When you select Any Legacy, the Legacy Data Converter examines all possible method templates in the source folder, regardless of version.

The conversion table displays the method templates in the Templates folder for the selected source version. The application verifies that the method template file is in the .pmtx file format.

3. To convert a method template that is not in the default list, do the following:

- a. Click **Browse** and locate a template folder.

You can select a specific template folder or a folder that contains multiple templates.

- b. Click **OK** in the Browse for Folder dialog box.

The application displays the selected folder in the conversion table.

When you select a folder that contains multiple method template folders, the application displays all the method templates.

4. In the Target Version list, select the version that you are converting to.

The list displays only TraceFinder configurations with compatible data. See [Version Compatibility](#).

5. (Optional) In the Target Name column, change the default name for each method template that you want converted.

When you populate the conversion table, the application checks each method template to see if a method template with this name exists in the target repository.

- If the method template name already exists in the target repository, the default new name appends “_1” to the original name.
- If the method template name does not exist in the target repository, the application keeps the original method template name.

IMPORTANT The conversion cannot overwrite an existing file name. If the new name is identical to an existing method template file, the conversion will fail. When you manually enter a new name, you must verify that the name does not already exist.

6. Select the **Convert** check box for each method template that you will convert, and click



The application confirms that all method templates to be converted use the .pmtx file format.

When the conversion process begins, the application displays a status bar and a Cancel button. You can cancel pending conversions, but not the template that is currently converting.

When the Status column reports that the template is successfully converted, the application writes the converted template to the specified target repository.

Note If a template conversion fails, the Status column displays “Conversion failed.” The log file contains details about the failed conversion.

7. To view a log of the conversion, click **View Log**.

The application opens a cumulative log file for the session in a Notepad text editor window.

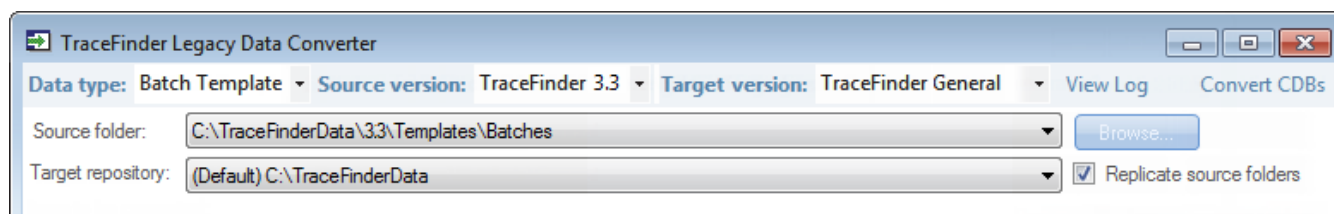
Converting Batch Templates

Use the data converter to convert legacy batch templates to TraceFinder 4.1 batch templates.

❖ To convert a batch template

1. In the Data Type list, select **Batch Template**.

The following example shows that you can convert batch templates from the TraceFinder 3.3 General configuration to the current General configuration. For a complete list of version compatibilities, see [Version Compatibility](#).



2. In the Source Version list, select the version of the batch template that you will convert.

Note When you select Any Legacy, the Legacy Data Converter examines all possible batch templates in the source folder, regardless of version.

The conversion table displays the batch templates in the Templates folder for the selected source version.

3. To convert a batch template that is not in the default list, do the following:
 - a. Click **Browse** and locate a template folder.

You can select a specific batch template folder or a folder that contains multiple batch templates.
 - b. Click **OK** in the Browse for Folder dialog box.

The application displays the selected folder in the conversion table.

When you select a folder that contains multiple batch template folders, the application displays all the batch templates.

4. In the Target Version list, select the version that you are converting to.

The list displays only TraceFinder configurations with compatible data. See [Version Compatibility](#).

5. (Optional) In the New Name column, change the default new name for each batch template that you want converted.

When you populate the conversion table, the application checks each batch template to see if a batch template with this name exists in the target repository.

- If the batch template name already exists in the target repository, the default new name appends “_1” to the original name.

- If the batch template name does not exist in the target repository, the application keeps the original batch template name.

IMPORTANT The conversion cannot overwrite an existing file name. If the new name is identical to an existing batch template file, the conversion will fail. When you manually enter a new name, you must verify that the name does not already exist.

6. Select the **Convert** check box for each batch template that you will convert, and click



The application confirms that all batch templates to be converted use the .btx file format.

When the conversion process begins, the application displays a status bar and a Cancel button. You can cancel pending conversions, but not the template that is currently converting.

When the Status column reports that the template is successfully converted, the application writes the converted template folder to the ...\\TraceFinderData\\4.0\\Templates\\Batches folder.

Note If a template conversion fails, the Status column displays “Conversion failed.” The log file contains details about the failed conversion.

7. To view a log of the conversion, click **View Log**.

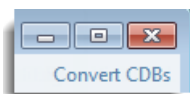
The application opens a cumulative log file for the session in a Notepad text editor window.

Converting Compound Databases

Use the data converter to convert legacy compound databases to TraceFinder 4.1 compound databases.

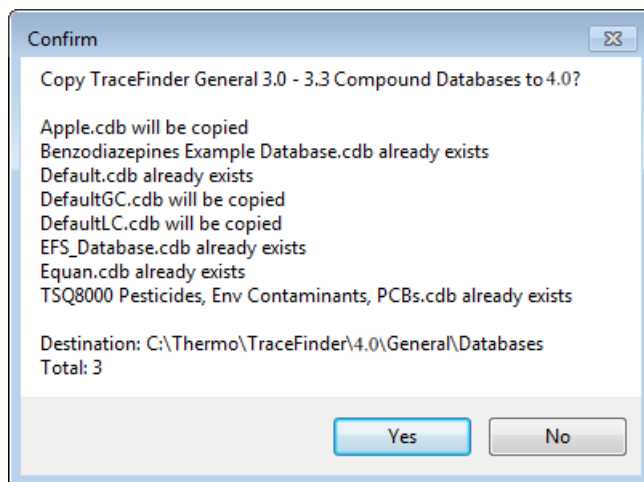
❖ To convert compound databases

1. Click **Convert CDBs** (in the upper right corner of the window).



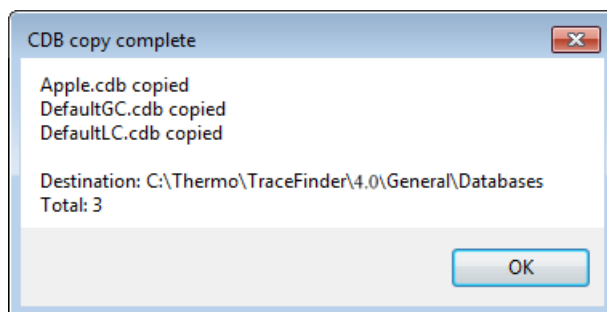
Note If your window is too narrow, you might not see the Convert CDBs button. Widen the window to expose the button in the upper right corner.

The Confirm dialog box lists all databases that will be converted to TraceFinder 4.1.



2. Click **Yes**.

The Confirm dialog box lists all databases that are converted to TraceFinder 4.1.



3. Click **OK**.

The converted compound databases are now available in the TraceFinder 4.1 application.

Choosing a Mode or Console

When user security is activated, the navigation pane displays the modes and consoles available to the current user's assigned roles and permissions. The following table shows the available modes and consoles for each user role.

Table 4. User roles and default access

User role	Method Development	Acquisition	Analysis	Configuration console	Administrator Console
Security					Security only
LabDirector	✓	✓	✓	✓	✓
ITAdmin					✓
Supervisor	✓	✓	✓	✓	✓
Technician		✓	✓		
QAQC			✓		

Note When user security is not activated, all modes and consoles are available to all users.

Follow these procedures:

- [To choose a mode](#)
- [To open the Configuration console](#)
- [To open the Administrator console](#)
- [To display a log of instrument errors](#)
- [To monitor instrument status](#)
- [To watch acquisition and processing in real time](#)

❖ **To choose a mode**


In the navigation pane, click the mode where you want to work.

The navigation pane shows only the modes that you have permission to use.



Mode	Description
Acquisition	Opens the Acquisition mode where you can create and review batches, batch data, reports, and local methods. See Chapter 3, “Using the Acquisition Mode.”
Analysis	Opens the Analysis mode where you can review batches, batch data, reports, and local methods. See Chapter 4, “Using the Analysis Mode for Quantitation Batches.” See Chapter 5, “Using the Analysis Mode for Target Screening Batches.” See Chapter 6, “Using the Analysis Mode for Unknown Screening Batches.”
Method Development	Opens the Method Development mode where you can create a master method or an instrument method. Refer to the <i>TraceFinder Lab Director User Guide</i> .

❖ **To open the Configuration console**

Click the **Application Configuration** icon, , in the upper right corner of the TraceFinder window.

When user security is activated, you must have Configuration permissions to access the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.

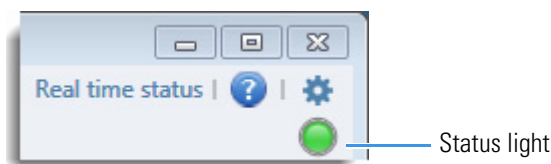
❖ **To open the Administrator console**

Choose **Tools > Administrator Console** from the TraceFinder main menu.

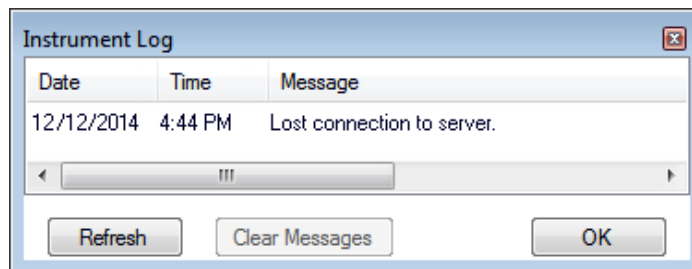
When user security is activated, you must have Administrator permissions to access the Administrator Console. Refer to the *TraceFinder Administrator Console User Guide*.

❖ To display a log of instrument errors

1. Click the status light in the upper right corner of the TraceFinder window.



The Instrument Log dialog box opens.



The Instrument Log displays all instrument errors that have occurred since the application started or since the last time that you cleared the message log.

2. Do any of the following:
 - Click **Refresh** to display errors that occur after you open the Instrument Log dialog box.
 - Click **Clear Messages** to remove messages from the Instrument Log display.

The application clears messages only from the Instrument Log display. These messages remain in the following log file:

C:\Thermo\TraceFinder\4.0\Clinical\Logs\TraceFinder.log

- Click **OK** to dismiss the Instrument Log dialog box.

❖ To monitor instrument status

Look at the status light in the upper right corner of the TraceFinder window.



Green indicates that the instrument is ready.





Yellow indicates that the instrument is in standby mode, waiting for contact closure, or warming up the lamp.



Red indicates that the instrument is not ready because it is turned off, not configured or connected, currently initializing, or otherwise busy.

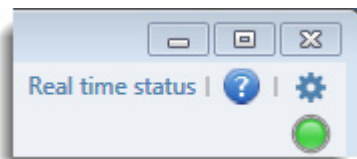
The status light (in any color) might include an exclamation mark.

- An exclamation mark on a green status light, , might indicate that the instrument method contains the wrong source (for example, APCI instead of HESI).
- An exclamation mark on a red status light, , might indicate that a vial is missing and the batch was forced to stop.

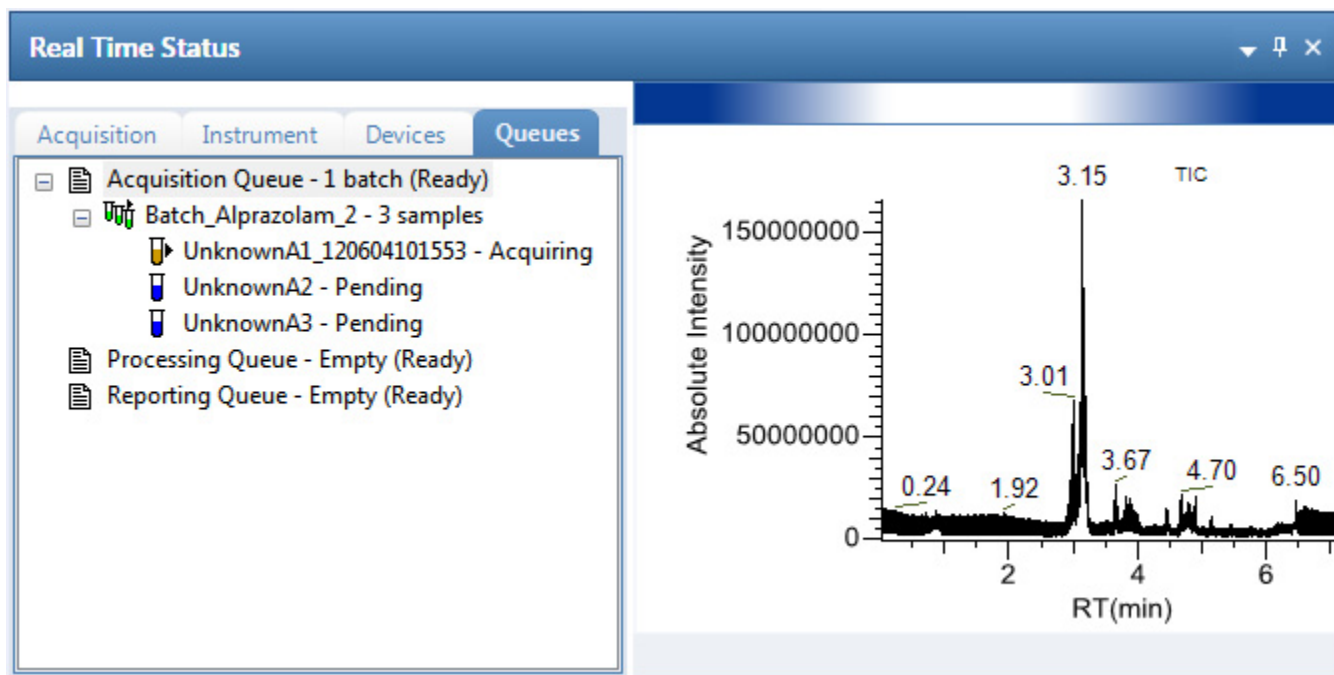
Click the status light to display details of the error.

❖ **To watch acquisition and processing in real time**

Click **Real Time Status** in the upper right corner of the TraceFinder window.



The application displays the Real Time Status pane at the bottom of the window.



For descriptions of all the features of the Real Time Status pane, see [Real Time Status Pane](#) in Chapter 3, "Using the Acquisition Mode."

Using the Acquisition Mode

This chapter describes the tasks associated with the Acquisition mode.

Contents

- [Working with Batches](#)
- [Real Time Status Pane](#)
- [Sample Types](#)

When you plan to work with multiple samples or use similarly designed batches, use the Acquisition mode to reduce the amount of data you must enter.

Because the nature and types of batches are often similar (in some cases specified by laboratory standard practices), you can define a batch template that supplies the basic structure of a batch.

IMPORTANT When user security is activated, you must have Acquisition Wizard – Template Editing permission to create a batch template.

Using a master method, you can create a batch and run the samples. A batch represents one or more samples that are to be acquired, processed, reviewed, and reported as a set. After you create a batch of samples, you can submit the batch and review the results in Data Review or you can go directly to viewing and printing reports.

You can set up a calibration batch with known concentrations of the target compounds and compare the calibration values against samples in future batches.

You can also use the Quick Acquisition feature to quickly submit samples from any page in the Acquisition mode. See [Appendix A, “Using Quick Acquisition.”](#)

Working with Batches

See the instructions for the following tasks:

- [Opening and Navigating the Acquisition Mode](#)
- [Creating and Submitting Batches](#)

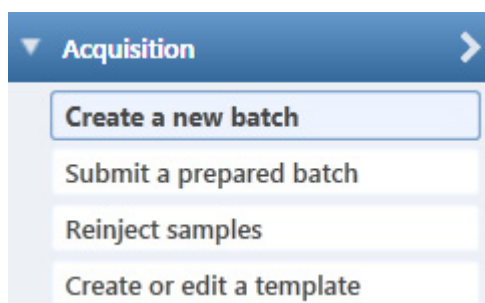
IMPORTANT TraceFinder 4.1 uses the same data as TraceFinder 4.0. By default, the application stores the batch data for the 4.1 release in the TraceFinderData\4.0\Projects folder.

Opening and Navigating the Acquisition Mode

❖ To access the Acquisition mode

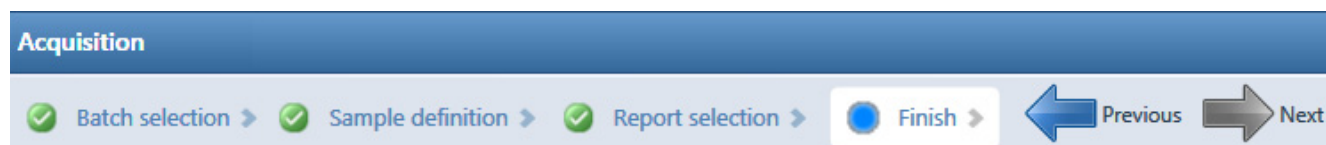
Click **Acquisition** in the navigation pane.

The navigation pane for the Acquisition mode opens.

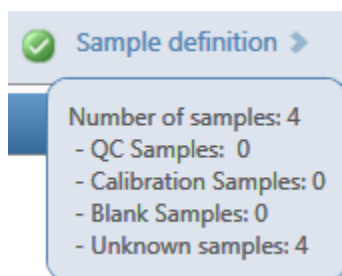
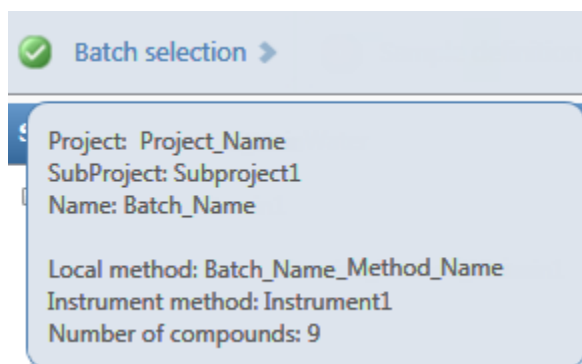


As you progress through the Acquisition mode using any of these methods for creating a batch, the task pane at the top of the view tracks your progress. As you complete each stage, you can hold your cursor over the view name in the task pane to display the parameters that you specified for the batch. See [Example task pane when you have completed the Acquisition mode](#).

Figure 6. Example task pane when you have completed the Acquisition mode



Hold your cursor over Batch Selection, Sample Definition, or Report Selection to view the parameters for your batch.



Categories in the Sample Definition list:

QC Samples: QC

Calibration Samples: Calibrator

Blank Samples: Negative

Unknown Samples: All other samples



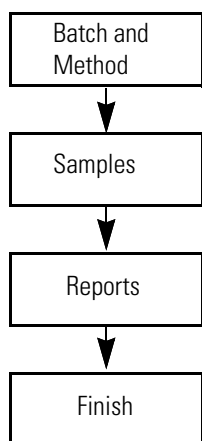
Creating and Submitting Batches

To create and submit a batch, the Acquisition mode uses a wizard-style interface to guide you through these major steps:

1. [Selecting a Batch](#)
2. [Defining a Sample List](#)
3. [Selecting and Reviewing Reports](#)
4. [Submitting a Batch in the Acquisition Wizard](#)

The Acquisition mode provides multiple techniques for creating either a batch or a batch template. Each batch creation technique has an associated workflow, as shown in the following flowcharts. Each workflow uses a different combination of Acquisition mode pages.

Workflow for Creating an Original Batch

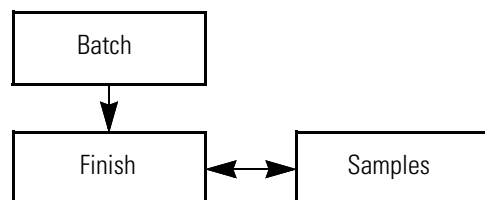


To create an original batch, start with any of these instructions: [To start a new quantitation batch](#), or [To start a new target screening batch](#), or [To start a new unknown screening batch](#).

❖ To see how to acquire a new batch

1. Choose **Help > Animations**.
2. From the list of animation topics, click **Acquiring a New Batch**.

Workflow for Acquiring a Prepared Batch

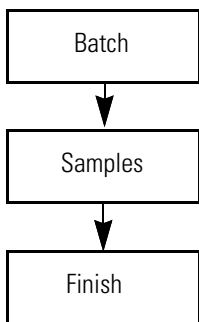


To acquire a prepared batch, start with the instructions [To select a prepared batch](#).

❖ **To see how to acquire a prepared batch**

1. Choose **Help > Animations**.
2. From the list of animation topics, click **Acquiring a Prepared Batch**.

Workflow for Reinjecting a Previously Acquired Batch

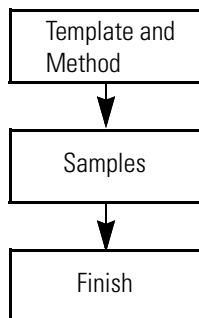


To process a previously acquired batch, start with the instructions [To reinject samples in a previously acquired batch](#).

❖ **To see how to reinject samples in a previously acquired batch**

1. Choose **Help > Animations**.
2. From the list of animation topics, click **Reinjecting Samples**.

Workflow for Creating or Editing a Batch Template



IMPORTANT When user security is activated, you must have Acquisition Wizard – Template Editing permission to create a batch template.

To create a batch template, start with any of these instructions: [To create a quantitation batch template](#), [To create a target screening batch template](#), or [To create an unknown screening batch template](#).

❖ **To see how to create a batch template**

1. Choose **Help > Animations**.
2. From the list of animation topics, click **Creating a Batch Template**.

Selecting a Batch

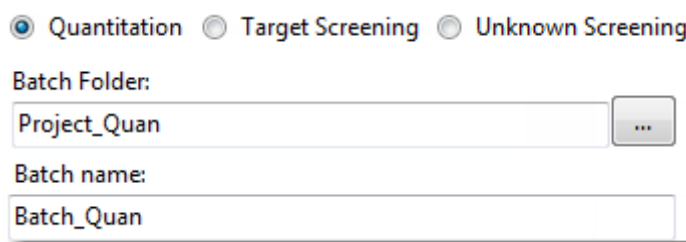
On the Batch Selection page of the Acquisition mode, you can create a new quantitative, target screening, or unknown screening batch in any of your current projects/subprojects. Or, you can submit a batch that you previously prepared and saved, reinject the samples in a batch that you previously acquired, or create a batch template to use for future batches.

Follow these procedures:

- [To start a new quantitation batch](#)
- [To start a new target screening batch](#)
- [To start a new unknown screening batch](#)
- [To start a new batch from a template](#)
- [To select a prepared batch](#)
- [To reinject samples in a previously acquired batch](#)
- [To create a quantitation batch template](#)
- [To create a target screening batch template](#)
- [To create an unknown screening batch template](#)

❖ To start a new quantitation batch

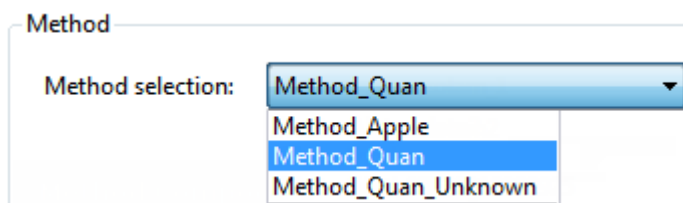
1. Click **Create a New Batch** in the navigation pane.
2. Select the **Quantitation** option.
3. Select the batch folder where you want to create the new batch.
4. Type a unique name for the new batch in the Batch Name box.



The screenshot shows a dialog box with three radio buttons at the top: **Quantitation** (selected), **Target Screening**, and **Unknown Screening**. Below the radio buttons, there are two text input fields. The first field is labeled "Batch Folder:" and contains the text "Project_Quan". To the right of this field is a button with three dots "...". The second field is labeled "Batch name:" and contains the text "Batch_Quan".

If the name you enter is not unique, a red warning flashes.

5. Select a method from the Method Selection list.



The screenshot shows a dropdown menu titled "Method". The label "Method selection:" is to the left of the dropdown. The dropdown is open, showing a list of methods: "Method_Quan" (highlighted in blue), "Method_Apple", "Method_Quan", and "Method_Quan_Unknown".

The Method Compound Data pane displays the compounds in the method. You cannot edit the compounds list from the Acquisition mode.

Method

Method selection: Method_Quan

Method Compound Data

	RT	Compound	Compound type	Active	CAS No
▶	3.14	Propanenitrile	Target Compound	<input checked="" type="checkbox"/>	107120
	3.15	EDIFENPHOS-CE5-R20-T...	Target Compound	<input checked="" type="checkbox"/>	17109498

- To continue to the next page, click **Next**.

The Sample Definition page opens. See [Defining a Sample List](#).

❖ To start a new target screening batch

- Click **Create a New Batch** in the navigation pane.
- Select the **Target Screening** option.
- Select the batch folder where you want to create the new batch.
- Type a name for the new batch in the Batch Name box.

☐ Quantitation
 ☒ Target Screening
 ☐ Unknown Screening

Batch Folder:

Project_Target ...

Batch name:

Batch_Target

If the name you enter is not unique, a red warning flashes.

- Select a method from the Method Selection list.

The Method Compound Databases pane displays the compound databases in the method. The application uses these databases to identify the compounds in the samples. You cannot edit the compound database list from the Acquisition mode.

Method

Method selection: Method_Screening

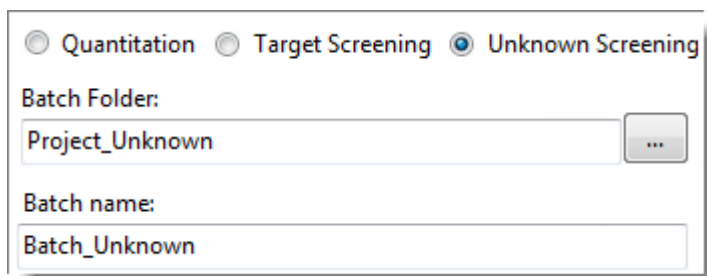
Method Compound Databases	
Enabled	Database File
▶ <input checked="" type="checkbox"/>	Default Database
<input checked="" type="checkbox"/>	General_Database
<input checked="" type="checkbox"/>	General_Database2

- To continue to the next page, click **Next**.

The Sample Definition page opens. See [Defining a Sample List](#).

❖ **To start a new unknown screening batch**

1. Click **Create a New Batch** in the navigation pane.
2. Select the **Unknown Screening** option.
3. Select the batch folder where you want to create the new batch.
4. Type a name for the new batch in the Batch Name box.



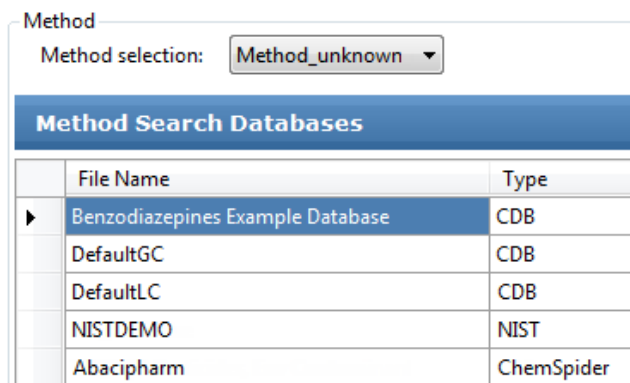
☐ Quantitation ☐ Target Screening ☒ Unknown Screening

Batch Folder:
Project_Unknown

Batch name:
Batch_Unknown

5. Select a method from the Method Selection list.

The Method Search Databases pane displays the databases that you selected in the method. The application uses these databases to identify the compounds in the samples. You cannot edit the database list from the Acquisition mode.



Method
Method selection: Method_unknown

Method Search Databases	
File Name	Type
▶ Benzodiazepines Example Database	CDB
DefaultGC	CDB
DefaultLC	CDB
NISTDEMO	NIST
Abacipharm	ChemSpider

6. To continue to the next page, click **Next**.

The Sample Definition page opens. See [Defining a Sample List](#).

❖ **To start a new batch from a template**

1. Click **Create a New Batch** in the navigation pane.
2. Select either the **Quantitation**, **Target Screening**, or **Unknown Screening** option.

The Available Templates pane displays only batch templates for the selected option.


3. In the Available Templates pane, select the template and method combination that you want to use.

Batch name:
Template_Quan_1

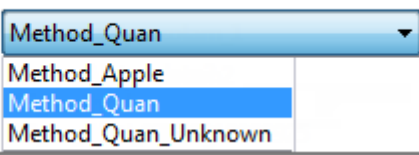
Available Templates

Template name	Method
Template_Quan	Method_Quan

The system creates a batch name with the selected template name and appends the date and time stamp. You can change the batch name, the default batch folder, or the method associated with this template.

4. (Optional) Click  and select a different batch folder where you want to create the new batch.
5. (Optional) Select a different method to use for the new batch.

Method

Method selection: 

Method_Quan
Method_Apple
Method_Quan
Method_Quan_Unknown

6. To continue to the next page, click **Next**.




The Sample Definition page of the Acquisition mode opens. See [Defining a Sample List](#).

❖ To select a prepared batch

1. Click **Submit a Prepared Batch** in the navigation pane.

The application displays all your unacquired, saved batches. The application stores all unacquired batches in the ...\\TraceFinderData\\4.0\\Projects\\... folder.

2. Select the batch that you want to acquire.

 Batch selection >  Previous  Next

Batch name	Folder	Last modified
Batch_Quan_1	Project_Quan	02-05-2015 11:00AM
Batch_Quan_1	Project_Quan	02-05-2015 10:59AM
Batch_Target_1	Project_Target	02-05-2015 11:00AM
Batch_Unknown_1	Project_Unknown	02-05-2015 11:01AM


3. To continue to the next page, click **Next**.

The Finish page of the Acquisition mode opens. From the Finish page, you can save the batch, submit the batch for acquisition, or go to the Sample Definition page to edit the sample list for this batch.

- If the batch is unreadable, the application reports that the batch file is not valid and cannot be opened.
 - If a sample in the batch is unreadable, the application cannot open the sample. The application creates a new sample with the same name and flags the sample. You must complete the missing information such as Sample Type, Level, and so forth, and then save the batch before you submit it for acquisition. Or, you can browse in a new raw data file to replace the corrupt file.
4. Do one of the following:
 - To edit the sample list, click **Previous**.


For detailed instructions, see [Defining a Sample List](#).

—or—

- To prepare the batch for acquisition, click **Submit**,  **Submit**.

For detailed instructions, see [Submitting a Batch in the Acquisition Wizard](#).

—or—

- To save the batch to be acquired later, click **Save**,  **Save**.

The application saves your batch in the following folder:

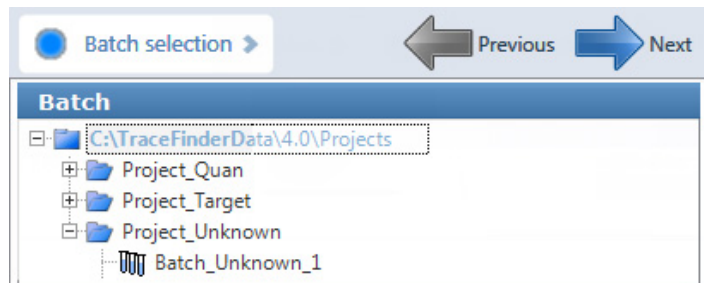
...\TraceFinderData\4.0\Projects\...

The application closes the Acquisition mode and returns you to the mode you were last using.

❖ **To reinject samples in a previously acquired batch**

1. Click **Reinject Samples** in the navigation pane.
2. On the Batch page, select the batch that you want to reacquire.

The Batch page displays all previously acquired quantitation, target screening, and unknown screening batches.



- To continue to the next page, click **Next**.

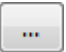
The Sample Definition page of the Acquisition mode opens. See [Defining a Sample List](#).

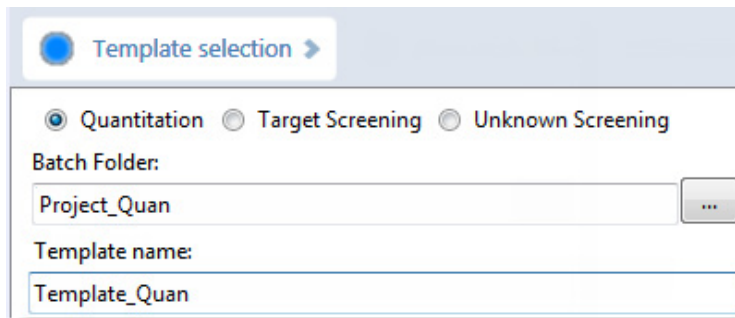
- If the batch is unreadable, the application reports that the batch file is not valid and cannot be opened.
- If a sample in the batch is unreadable, the application creates a new sample with the same name and flags the sample. You must complete the missing information, such as Sample Type, Level, and so forth, and then save the batch before you submit it for acquisition. Or, you can browse in a new raw data file to replace the corrupt file.

❖ To create a quantitation batch template

- Click **Create or Edit a Template** in the navigation pane.

Note When user security is activated, you must have Acquisition Wizard – Template Editing permission to create a batch template.

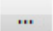
- Select the **Quantitation** option.
- Click  and select the folder where you want to create the new batch template.
- Type a name for the new batch template in the Template Name box.



Template selection >

☒ Quantitation ☐ Target Screening ☐ Unknown Screening

Batch Folder:

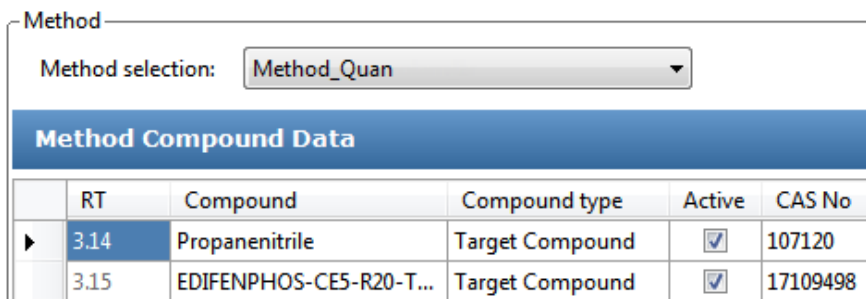
Project_Quan 

Template name:

Template_Quan

- Select a method from the Method Selection list.

The Method Compound Data pane displays the compounds in the method. You cannot edit the compounds list from the Acquisition mode.



Method

Method selection: Method_Quan

Method Compound Data					
	RT	Compound	Compound type	Active	CAS No
▶	3.14	Propanenitrile	Target Compound	<input checked="" type="checkbox"/>	107120
	3.15	EDIFENPHOS-CE5-R20-T...	Target Compound	<input checked="" type="checkbox"/>	17109498

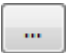
- To continue to the next page, click **Next**.

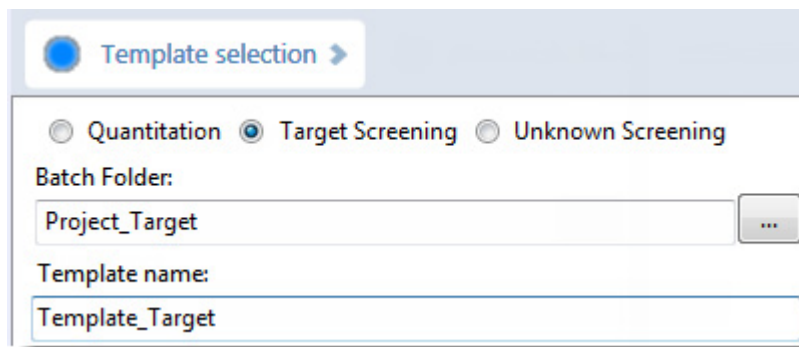
The Sample Definition page of the Acquisition mode opens. See [Defining a Sample List](#).

❖ **To create a target screening batch template**

1. Click **Create or Edit a Template** in the navigation pane.

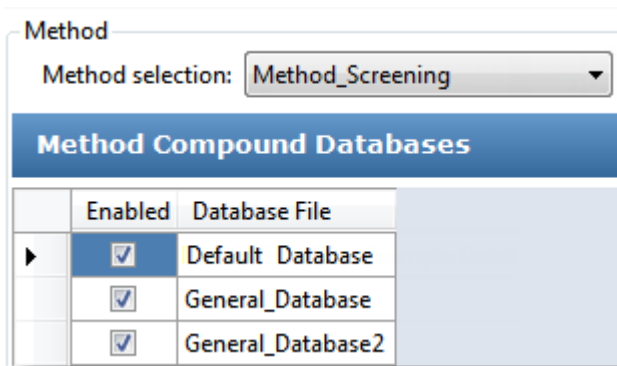
Note When user security is activated, you must have Acquisition Wizard – Template Editing permission to create a batch template.

2. Select the **Target Screening** option.
3. Click  and select the folder where you want to create the new batch template.
4. Type a name for the new batch template in the Template Name box.



5. Select a method from the Method Selection list.

The Method Compound Databases pane displays the screening databases available for the selected method.



	Enabled	Database File
▶	<input checked="" type="checkbox"/>	Default Database
	<input checked="" type="checkbox"/>	General_Database
	<input checked="" type="checkbox"/>	General_Database2

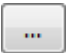
6. Select the check box for each database that you want to use for screening.
7. To continue to the next page, click **Next**.

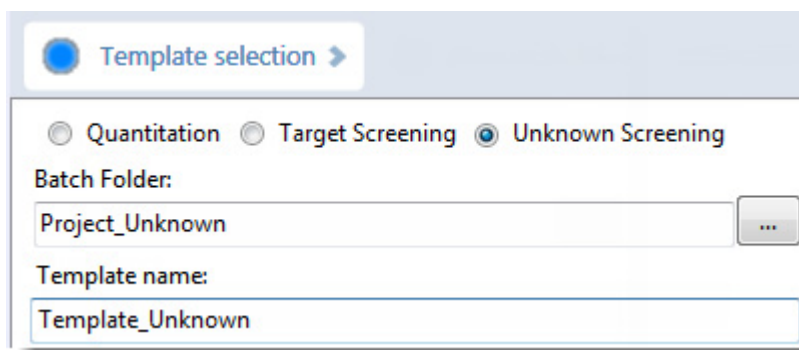
The Sample Definition page of the Acquisition mode opens. See [Defining a Sample List](#).

❖ **To create an unknown screening batch template**

1. Click **Create or Edit a Template** in the navigation pane.

Note When user security is activated, you must have Acquisition Wizard – Template Editing permission to create a batch template.

2. Select the **Unknown Screening** option.
3. Click  and select the folder where you want to create the new batch template.
4. Type a name for the new batch template in the Template Name box.



5. Select a method from the Method Selection list.

The Method Compound Databases pane displays the unknown screening databases available for the selected method.

6. Select the check box for each database that you want to use for unknown screening.
7. To continue to the next page, click **Next**.

The Sample Definition page of the Acquisition mode opens. See [Defining a Sample List](#).

Defining a Sample List

Use the [Samples](#) page on the Sample Definition page of the Acquisition mode to create a list of samples for the batch. You can add samples, insert samples, import a sample list, or remove samples from the list. You can use the [Reference Sample](#) page to select a reference sample to use as a reference peak in the Data Review.

To create the sample list, you can use either of two sets of function buttons (described in the following graphic) or you can use commands in the shortcut menu (see the Shortcut menu commands area of the [Samples page parameters](#)).

As you enter sample values, you can use the Copy Down and Fill Down commands to quickly enter column values. For detailed instructions on using Copy Down and Fill Down to enter column values, see [Appendix B, “Using Copy Down and Fill Down.”](#)




Use any of the following procedures to create a sample list. When you finish defining the list of samples, click **Next**.

- When you create a batch from scratch and click Next, the Report Selection page opens. See [Selecting and Reviewing Reports](#).
- When you edit a prepared batch, reinject samples, create a batch from a template, or edit a batch template and then click Next, the Finish Selection page opens. See [Submitting a Batch in the Acquisition Wizard](#).

Follow these procedures:

- [To add samples to the list](#)
- [To insert samples into the list](#)
- [To import samples into the list](#)
- [To remove samples from the list](#)
- [To reinject a sample from a previously acquired batch](#)
- [To select channels for the batch](#)
- [To assign a specific channel to a sample](#)
- [To select a reference sample](#)
- [To add an auto sample type](#)
- [To specify different instrument methods for samples](#)

❖ **To add samples to the list**

1. Select the number of sample rows to add  and then click the **Add** icon,  or .
2. Type a file name in the Filename column for each sample.
Each file name must be unique.
3. Select a sample type from the Sample Type list for each sample.

Available sample types

Specimen	Hydrolysis	Solvent	QC
Unextracted	Calibrator	Negative	

For a detailed description of each sample type, see [Sample Types](#).

4. For each Calibrator or QC sample, select a level from the Level list.
The master method defines the sample levels. If there are no levels to select in the Level list, ask a user with Method Development permission to edit the method and specify the levels. Then return to the Acquisition mode, and begin the batch again. The application does not save a batch when you leave the Acquisition mode.




If you have Method Development permission, do the following:

- a. Return to the Method Development mode.
- b. Open the method.
- c. Click the **Compounds** tab.
- d. Click the **Calibration Levels** tab.
- e. Add the levels.
- f. Save the method.

For detailed instructions, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

5. (Optional) To assign samples to a group, do the following:
 - a. For each sample, type a group name in the Groups column.
Repeat this for each sample that you want to include in a group.
 - b. Create as many groups as you want.

Note To assign a sample to multiple groups, separate the groups with a comma.

Status	Filename	Groups
	Benzo26473	groupB, groupA
	Benzo25557	groupB
	Benzo26154	groupB, groupA

The application uses only the first group listed for calculating the Maximum Fold and Group Averages.

For information about Maximum Fold analysis in unknown screening analysis, see [Cross Sample Peak List Pane](#).

- For quantitation experiments, after creating groups, you can choose one of the samples as a threshold sample for the group and then compare all of the samples in the Comparative View in Data Review.

For information about specifying a threshold sample for a group of samples, see [Threshold Samples Page for Quantitation Batches](#). For information about viewing grouped samples in Data Review, see [Comparative View for Quantitation Batches](#). For information about Group Averages, see [Group Averages](#).

- For unknown screening experiments, you can create a control group (specify the group name as Control) to use for group averages in Data Review. For information about Group Averages, see [Group Averages Pane](#).

- For each sample, type a vial position in the Vial Position column.

Tip Use the Fill Down command to make entering vial positions easier.

- For each sample, type a volume in the Injection Volume column.

The minimum injection volume value allowed is 0.1 µL; the maximum injection volume value allowed is 5000 µL.

- (Optional) Type or edit the values for the remaining columns.


Note When you use the scroll bar at the bottom of the sample list, the following columns remain fixed: Status, Filename, Sample Type, Groups, Qual Processing (quantitation) or Blank Subtraction (target screening), Level, Sample ID, and Sample Name, while the other columns scroll right and left.

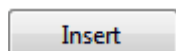
To automatically copy or fill values in these columns, see [Appendix B, “Using Copy Down and Fill Down.”](#)

❖ **To insert samples into the list**





1. Select the sample above which you want to insert new Specimen samples.

You cannot use the Insert command to create the first sample row.

2. Select the number of samples to insert and then click the **Insert** icon,  or



The application inserts the Specimen samples above the selected sample.

	Status	Filename	Sample type	Groups
1		cal_std_5	Calibrator	
2		Unknown2	Specimen	
3		Unknown1	Specimen	
4		cal_std_10	Calibrator	

3. For each sample, type a file name in the Filename column.
Each file name must be unique.
4. For each sample, select a sample type from the Sample Type list.

Available sample types

Specimen	Hydrolysis	Solvent	QC
Unextracted	Calibrator	Negative	




5. For each Calibrator or QC sample, click the Level cell and select a level from the list.

The master method defines the sample levels. If there are no levels to select from the Level list, ask a user with Method Development permission to edit the method and specify the levels. Then return to the Acquisition mode, and begin the batch again. The application does not save a batch when you leave the Acquisition mode.

If you have Method Development permission, follow the instructions in [step 4](#) of the procedure [To add samples to the list](#).

6. (Optional) To assign samples to a group, do the following:
 - a. For each sample, type the group name in the Groups column.
Repeat this for each sample that you want to include in a group.
 - b. Create as many groups as you want.

Note To assign a sample to multiple groups, separate the groups with a comma.

Status	Filename	Groups
	Benzo26473	groupB, groupA
	Benzo25557	groupB
	Benzo26154	groupB, groupA

The application uses only the first group listed for calculating the Maximum Fold and Group Averages.

For information about Maximum Fold analysis in unknown screening analysis, see [Cross Sample Peak List Pane](#).

- For quantitation experiments, after creating groups, you can choose one of the samples as a threshold sample for the group and then compare all of the samples in the Comparative View in the Analysis mode.

For information about specifying a threshold sample for a group of samples, see [Threshold Samples Page for Quantitation Batches](#). For information about viewing grouped samples in Data Review, see [Comparative View for Quantitation Batches](#). For information about Group Averages, see [Group Averages](#).

- For unknown screening experiments, you can create a control group (specify the group name as Control) to use for group averages in Data Review. For information about group averages, see [Group Averages Pane](#).

- Type a vial position in the Vial Position column for each sample.

Tip Use the Fill Down command to make entering vial positions easier.

- For each sample, type a volume in the Injection Volume column.

The minimum injection volume allowed is 0.1 µL; the maximum injection volume allowed is 5000 µL.

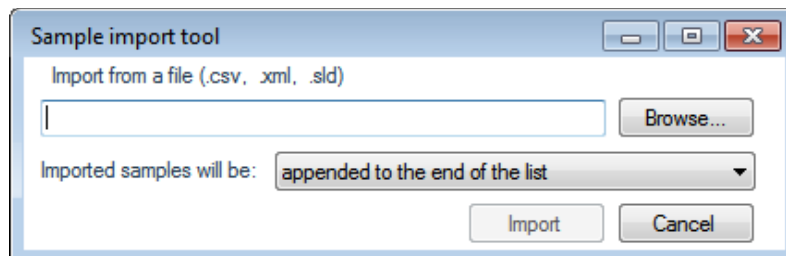
- (Optional) Type or edit the values for the remaining columns.

Note When you use the scroll bar at the bottom of the sample list, the following columns remain fixed: Status, Filename, Sample Type, Groups, Qual Processing (quantitation) or Blank Subtraction (target screening), Level, Sample ID, and Sample Name, while the other columns scroll right and left.

❖ **To import samples into the list**

1. Click **Import**, .

The Sample Import Tool dialog box opens.



Use this dialog box to import a sample list from a CSV, an XML, or an SLD file.

2. Click **Browse** and select a CSV, an XML, or an SLD file with the sample definitions that you want to import.

Note The .csv, .xml, or .sld data format must match the TraceFinder data format.

3. From the Imported Samples Will Be list, select either **Appended to the End of the List** or **Inserted at the Selected Row**.
4. Click **Import**.

The Sample Import Tool dialog box closes, and the application adds the specified samples to the sample list.

When you import samples from an Xcalibur sequence file (.sld), the application makes the following column name substitutions.

Xcalibur column	TraceFinder column
Position	Vial position
Inj Vol	Injection volume
Dil Factor	Conversion Factor

When you import samples from an Xcalibur sequence file (.sld), the application makes the following sample type substitutions.

Xcalibur sample type	TraceFinder sample type
Blank	Negative
Std Bracket	Calibrator

For each imported sample, the application uses the Instrument Method specified in the local method.

5. For each Calibrator or QC sample, click the Level cell and select a level from the list.

The master method defines the sample levels. If there are no levels to select from the Level list, ask a user with Method Development permission to edit the method and specify the levels. Then return to the Acquisition mode, and begin the batch again. The application does not save a batch when you leave the Acquisition mode.

If you have Method Development permission, follow the instructions in [step 4](#) of the procedure [To add samples to the list](#).

For detailed instructions about defining calibration levels, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

6. Type a vial position in the Vial Position column for each sample.

Tip Use the Fill Down command to make entering vial positions easier.

7. Type a volume in the Injection Volume column for each sample.

The minimum injection volume value allowed is 0.1 µL; the maximum injection volume value allowed is 5000 µL.

8. (Optional) Type or edit the values for the remaining columns.

Note When you use the scroll bar at the bottom of the sample list, the following columns remain fixed: Status, Filename, Sample Type, Groups, Qual Processing (quantitation) or Blank Subtraction (target screening), Level, Sample ID, and Sample Name, while the other columns scroll right and left.

9. (Optional) When using multiplexing, select a channel for each imported sample.
Imported samples default to Auto.

❖ **To remove samples from the list**

1. Select the samples that you want to remove.

Tip Use the CTRL or SHIFT keys to select multiple samples.

2. Right-click and choose **Remove Selected Samples**.





❖ **To reinject a sample from a previously acquired batch**

1. In the sample list, select the sample to reinject.
2. Right-click and choose **Reinject Selected Samples**.

The application creates a copy of the selected sample and appends INJ001 to the file name. Additional reinjections of the same sample are numbered INJ002, INJ003, and so forth.

The application copies all parameter values from the original sample.

A green status icon indicates previously acquired samples (acquired and processed), and the sample name is grayed out. A blue status icon indicates samples created for reinjection (not acquired).

	cal_50_INJ001	Calibrator	10
	cal_50	Calibrator	10
	cal_10_INJ001	Calibrator	10
	cal_10	Calibrator	10

When you submit this batch, the application acquires only the reinjection samples.

❖ To select channels for the batch

Note These features are available only when you have activated multiplexing in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.

To disable a configured channel, clear the check box for the channel in the Multiplexing Channels area at the bottom of the page.

Multiplexing Channels

☒ All Channels ☒ Channel 1 ☒ Channel 2 ☒ Channel 3 ☒ Channel 4

By default, all configured channels are selected. The configured channels are determined by the multiplexing settings in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.

Clearing a channel in the Multiplexing Channels area does not remove this channel selection from the Channels list for each sample. When you assign a channel to a sample, be careful not to assign a channel that is not available.

❖ To assign a specific channel to a sample

1. Scroll to the Channel column.

Note The Channel column is available only when you have activated multiplexing in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.

All samples default to Auto.

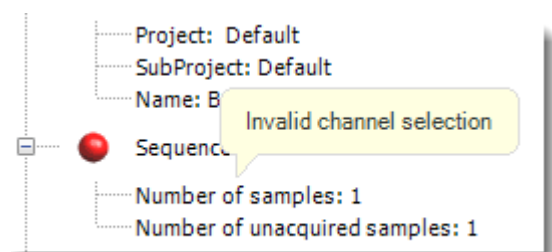
2. Select a channel from the Channel list.

Auto ▼

- Auto
- Channel 1
- Channel 2
- Channel 3
- Channel 4

When you submit the batch, samples that are set to Auto run on any of the available channels and samples that are set to a specific channel run only on that channel.

If you select a channel that is not available for this batch, the application flags the sample sequence on the Finish page of the Acquisition mode. See the previous procedure, [To select channels for the batch](#).



3. If you see this error, do the following:
 - a. Click **Previous** to return to the Sample Definition page.
The incorrect sample is marked with an error flag.
 - b. Correct the channel selection.

❖ To select a reference sample

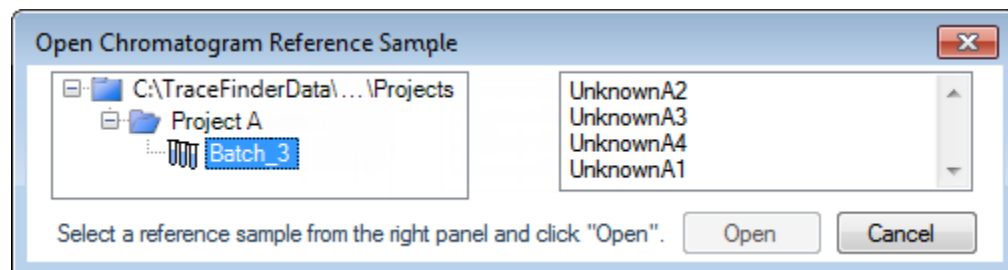
1. Click the **Reference Sample** tab.

The [Reference Sample](#) page on the Sample Definition page opens.

You can select one reference sample to use as a reference peak in the Data Review.

2. Right-click the Reference Sample page and choose **Add Reference Sample**.

The Open Chromatogram Reference Sample dialog box opens.



Note If you are using a new method, you will not see any samples here. You must create and save a batch using the current method to see available samples in this list.

3. Select a batch from the list.

The application displays only batches that were created using the current master method.

4. Select a sample from the list of processed samples on the right.

The application displays all the processed samples in the selected batch. To use a sample as a reference sample, the sample must have been processed with the current master method.

5. Click **Open**.

The application adds the reference sample to the Reference Sample page.

6. (Optional) Enter values for Sample ID, Sample Name, Comment, and Barcode Actual.
7. (Optional) Change the Vial Position for the sample.

The application uses the peak in this sample as a reference peak in Data Review. See [Reference Peak](#).

❖ To add an auto sample type

1. Click the **Auto Samples** tab.

The [Auto Samples](#) page opens.

2. Right-click and choose **Add Auto Sample**, or click the **Add New Auto Sample** icon,



The application adds a Solvent sample to the sample list.

You can add, insert, or remove samples from this list as you would any sample list.

3. To change the sample type to a Negative, click the Sample Type column and select Negative from the list.
4. In the Injection Volume column for the sample, type a volume.

The minimum injection volume value allowed is 0.1 µL; the maximum injection volume value allowed is 5000 µL.


5. In the Number of Injections column, type the number of injections available in the designated Solvent or Negative vial.

After auto sample injections have occurred, you can return to this page to view the number of Injections Used in each vial.

6. In the Vial Position column, type the vial position for the Solvent or Negative sample.

❖ **To specify different instrument methods for samples**

Note By default, the Instrument Method column is not displayed on the Sample Definition page. See [Instrument method column](#).

1. Display the Instrument Method column in the sample list:
 - a. Right-click the sample list and choose **Modify Columns**.
The Modify Columns dialog box opens.
 - b. In the Available Columns pane, select **Instrument Method**.
 - c. Click  to move the Instrument Method column to the Displayed Columns pane.
 - d. Click **OK**.

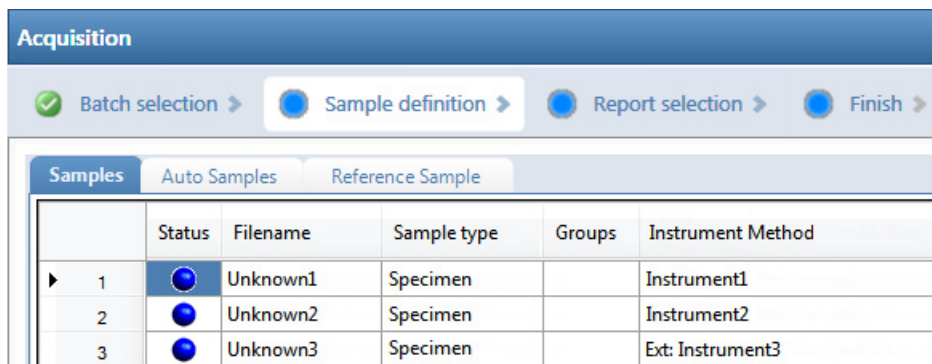
The application displays the Instrument Method column, defaulting to the instrument method specified in the master method.




2. Click the Instrument Method column and select an instrument method from the list.

This list contains all the available instrument methods. The application prefixes instrument methods from external sources with “Ext:”.

You can specify a different instrument method for each sample.

Figure 7. Instrument method column



	Status	Filename	Sample type	Groups	Instrument Method
1		Unknown1	Specimen		Instrument1
2		Unknown2	Specimen		Instrument2
3		Unknown3	Specimen		Ext: Instrument3

When you submit the batch, the application saves a copy of the selected instrument methods to the following folders:

External instrument methods:

...\\TraceFinderData\\4.0\\Projects\\...\\batch\\Methods\\method\\ExternalMethods

Local instrument methods:

...\\TraceFinderData\\4.0\\Projects\\...\\batch\\Methods\\method

Samples

Use the features on the Samples page to create a list of samples for the batch.

Figure 8. Samples page on the Sample Definition page

The screenshot shows the 'Samples' page within the 'Sample Definition' section. At the top, there are navigation buttons: 'Batch selection', 'Sample definition' (active), 'Report selection', and 'Finish'. Below these are 'Previous' and 'Next' buttons. The main area contains a table with the following data:

	Status	Filename	Sample type	Groups	Level	Sample ID	Sample name	Vial position	Injection volume	Channel	Barcode Expected	Sample Volume	Dilution Factor	Sample Weight	Calculation Type
1		Unknown1	Solvent					CStk1-01:1	10.0	Channel 1		1	1	1	Liquid
2		Unknown2	Negative					CStk1-01:3	10.0	Auto		1	1	1	Liquid
3		Unknown3	Specimen					CStk1-01:4	10.0	Auto		1	1	1	Liquid

Below the table, there are 'Sample Controls' with 'Add', 'Insert', and 'Import' buttons, each with a numeric input field. To the right, 'Multiplexing Channels' are listed with checkboxes: 'All Channels', 'Channel 1', 'Channel 2', 'Channel 3', and 'Channel 4'.

Table 5. Samples page parameters (Sheet 1 of 2)

Parameter	Definition
Previous	Returns you to the previous Acquisition page.
Next	Takes you to the next Acquisition page.
Status	Sample is not acquired. Sample is acquired but not processed. Sample is acquired and processed. Sample is currently acquiring.
Sample Controls	
Add	Adds the specified number of empty rows to the sample grid.
Insert	Inserts the specified number of empty rows above the selected row.
Import	Opens the Sample Import Tool to import samples from a CSV, an XML, or an SLD file.
Multiplexing Channels	These features are available only when you have activated multiplexing in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the <i>TraceFinder Lab Director User Guide</i> .
All Channels	Uses all configured channels to acquire this batch.
Channel 1- <i>n</i>	Uses only the selected channels to acquire this batch.
Shortcut menu commands	
Add Sample	Adds a single empty row to the sample grid.
Insert Sample	Inserts a single empty row to the sample grid above the selected row.
Insert Copy Sample	Copies the currently selected row and inserts a copy above the row.
Reinject Selected Samples	Creates a copy of the selected sample and appends INJ001 to the file name. Additional reinjections of the same sample are numbered INJ002, INJ003, and so forth.

Table 5. Samples page parameters (Sheet 2 of 2)

Parameter	Definition
Remove Selected Samples	Removes selected samples from the sample grid.
Import Samples	Opens the Sample Import Tool. Follow the instructions To import samples into the list.
Copy Down	Copies the value in the selected row to all rows below it. For detailed instructions about using the Copy Down command, see Appendix B, “Using Copy Down and Fill Down.”
Fill Down	Enters sequential values in the column starting with the value in the selected row and ending with the last row in the column. For detailed instructions about using the Fill Down command, see Appendix B, “Using Copy Down and Fill Down.”
Modify Columns	Opens the Modify Columns dialog box. See Column Display .
Enable/Disable Sample Weight Calculation	Displays or hides the Sample Volume, Dilution Factor, Sample Weight, Calculation Type, and Final Units columns.
Copy	Copies the data in the selected rows or columns to the Clipboard. Use this command to copy sample information to a text editor or spreadsheet application. You cannot paste this data back into the Acquisition mode sample list.
Copy with Headers	Copies the data in the selected rows or columns and the associated column headers to the Clipboard. Use this command to copy sample information to a text editor or spreadsheet application. You cannot paste this data back into the Acquisition mode sample list.
Paste	Pastes a single column of copied data from a text editor or spreadsheet application, into the selected column.
Undo Last Paste	Removes the last pasted item in the Acquisition mode sample list.
Export to CSV File	Opens the Save As dialog box where you can save the current sample list to a CSV file.
Edit Instrument Method	<p>Opens the Instrument Setup window where you can edit the parameters of the instrument method.</p> <ul style="list-style-type: none"> • When you edit an external method, the application updates the method in the ...\\Xcalibur\\methods folder. • When you edit an internal method, the application updates the method in the ...\\TraceFinderData\\4.0\\Projects\\<i>project\\subproject\\batch</i>\\Methods\\<i>method</i> folder. <p>For detailed information about editing instrument methods, refer to Chapter 3, “Using Instrument Methods in the Method Development Mode,” in the <i>TraceFinder Lab Director User Guide</i>.</p>

Column Display

The Samples page contains many columns of information. You can scroll to see all the columns, and you can customize which ones to display and their display order.

❖ To scroll the sample list

Use the horizontal scroll bar at the bottom of the sample list to view all the information.

When you use the scroll bar at the bottom of the sample list, the following columns remain fixed: Status, Filename, Sample Type, Groups, Qual Processing (quantitation) or Blank Subtraction (target screening) Level, Sample ID, and Sample Name, while the other columns scroll right and left.

❖ To customize the column display

1. Right-click the sample list and choose **Modify Columns**.

The [Modify Columns dialog box](#) opens.

2. Use the arrow buttons to move all the columns that you want displayed to the Displayed Columns pane.

These columns appear after the Status, Filename, Sample Type, Groups, Qual Processing (quantitation) or Blank Subtraction (target screening), Level, Sample ID, and Sample Name columns.

3. To arrange the order of the columns, do the following:
 - a. In the Displayed Columns pane, select a column name.
 - b. Use **Up** or **Down** to move the selected column up or down in the list.

The first column in the list represents the leftmost column in the Samples page sample list, and the last column in the list represents the rightmost column in the Samples page sample list.

Note The following columns are fixed: Status, Filename, Sample Type, Groups, Qual Processing (quantitation) or Blank Subtraction (target screening), Level, Sample ID, and Sample Name.

4. To change the width of a column, do the following:
 - a. In the Displayed Columns pane, select the column width.

5	Sample ID	100
▶ 6	Sample name	100
7	Vial position	100

- b. Type a new value for the width.
5. Repeat step 4 for all columns whose widths you want to change, and click **OK**.

The columns in the sample list immediately reflect your changes.

Figure 9. Modify Columns dialog box

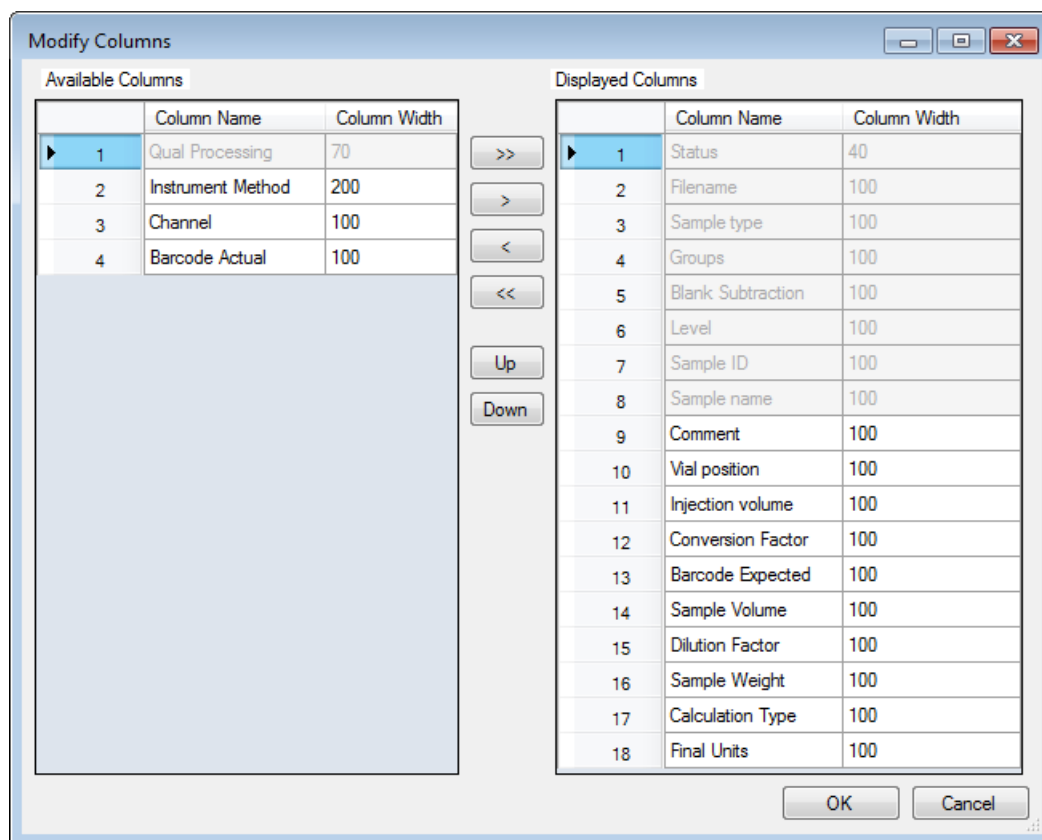



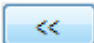
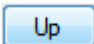
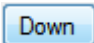


Table 6. Button descriptions for the Modify Columns dialog box

Button	Description
	Moves all columns to the Displayed Columns pane.
	Moves the selected column to the Displayed Columns pane.
The following buttons apply to all columns, except for those that are fixed: Status, Filename, Sample Type, Groups, Qual Processing (quantitation) or Blank Subtraction (target screening), Level, Sample ID, and Sample Name.	
	Moves the selected column to the Available Columns pane.
	Moves all columns except fixed columns.
	Moves the selected column name in the Displayed Columns pane one row up in the column order.
	Moves the selected column name in the Displayed Columns pane one row down in the column order.

Auto Samples

Use the features on the Auto Samples Sample page to identify the Solvent or Negative samples to use for any Auto Sample or Auto Sample and Reinject failure actions as specified on the Intelligent Sequencing page of the method. Refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

Each sample type that you specify for a failure action on the Intelligent Sequencing page must be defined in the samples list on the Auto Samples page.

Figure 10. Auto Samples page on the Sample Definition page

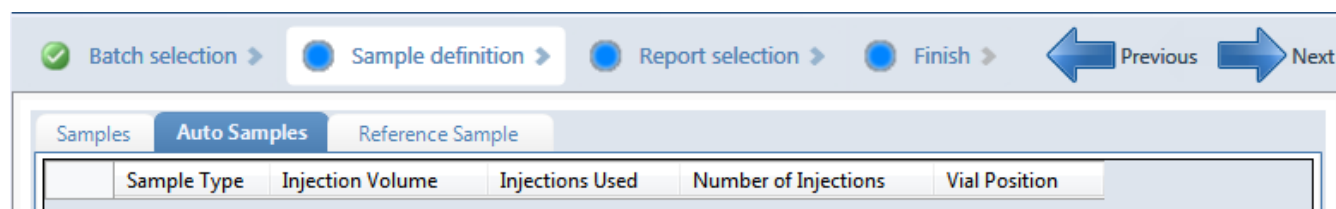


Table 7. Auto Samples page parameters

Column	Description
Sample Type	The sample type for the auto sample injection as specified on the Intelligent Sequencing page of the method—either Solvent or Negative. Default: Solvent
Injection Volume	The injection volume used for the sample acquisition as specified on the Samples page. Valid range: 0.1 through 5000 µL
Injections Used	The number of times a vial has been used. The count is cumulative across all batches.
Number of Injections	The number of injections available in the designated Solvent or Negative vial.
Vial Position	Vial position for this sample type as specified on the Samples page.

Reference Sample

Use the features on the Reference Sample page to select a sample to use as a reference peak in Data Review.

Figure 11. Reference Sample page on the Sample Definition page

	Status	Filename	Sample ID	Sample name	Comment	Vial position	Barcode Actual
▶ 1		Unknown1				1	

Table 8. Reference Sample page parameters

Parameter	Description
Status	Sample is not acquired. Sample is acquired but not processed. Sample is acquired and processed. Sample is currently acquiring.
Filename	Name of the raw data file that contains the sample data.
Sample ID	A user-defined, alphanumeric string that identifies a sample.
Sample Name	A user-defined name that identifies a sample.
Vial Position	The tray vial number used for an autosampler acquisition.
Barcode Actual	A user-entered barcode for the vial.
Shortcut menu commands	
Add Reference Sample	Opens the Open Chromatogram Reference Sample dialog box where you can select a reference sample.
Delete Selected	Deletes the reference sample.
Copy	Copies the data in the selected rows or columns to the Clipboard. Use this command to copy sample information to a text editor or spreadsheet application. You cannot paste this data back into the reference sample list.
Copy with Headers	Copies the data in the selected rows or columns and the associated column headers to the Clipboard. Use this command to copy sample information to a text editor or spreadsheet application. You cannot paste this data back into the reference sample list.
Paste	Pastes a single column of copied data from a text editor or spreadsheet application, into the selected column.
Export to CSV File	Opens the Save As dialog box where you can save the current sample list to a CSV file.

Selecting and Reviewing Reports

On the Report Selection page, you can specify the types of reports that you want to create. See [Report Selection](#). In addition to the report type, you can specify a report description for each of your reports.

For each report that you generate, you can create a hard-copy printout, a PDF file, a CSV file, or an Excel file.

Use any of the following procedures to create a reports list. When you finish specifying your report options, click **Next** to go to the Finish page and submit your batch. See [Submitting a Batch in the Acquisition Wizard](#).

The application writes the resulting output files for your reports to the following folder:

...\\TraceFinderData\\4.0\\Projects\\...*batch*\\Reports

Follow these procedures:

- [To edit a report title](#)
- [To specify a report in print format or as a PDF, a CSV, or an Excel file](#)

❖ To edit a report title

Select the Report Title column and edit the default title.

The default report title is the same as the report name.

❖ To specify a report in print format or as a PDF, a CSV, or an Excel file

1. For each type of report that you want to create, select the corresponding check box in the Print, Create PDF, Create CSV, or Create Excel column.
2. To duplicate the output type for all reports, right-click the cell and choose **Copy Down**.

All check boxes in the column below the selected cell duplicate the selected or cleared state in the selected cell. This action applies only to report types that make this output format available.

Report Selection

Use the features on the Report Selection page to specify the types of reports that you want to create.

Figure 12. Report Selection page

Report Name	Report Title	Print	Create PDF	Create CSV	Create Excel
Batch Report	Batch Report	<input type="checkbox"/>	<input type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>
Blank Report	Blank Report	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
Calibration Report	Calibration Report	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
Check Standard Report	Check Standard Report	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>

Table 9. Report Selection page parameters

Parameter	Description
Report Name	The name of a report.
Report Title	User-editable description to be used on a report.
Print	Reports to be sent to the printer.
Create PDF	Reports to be saved as PDF files.
Create CSV	Reports to be exported as CSV files.
Create Excel	Reports to be exported as Excel files.
Shortcut menu: Copy Down	Copies the selected or cleared state to all subsequent reports in the column.

Submitting a Batch in the Acquisition Wizard

On the [Finish](#) page of the Acquisition mode, you can specify a startup method, a shutdown method, or a calibration batch. You can save the batch to be acquired later, or you can acquire and process data and optionally create reports.

Note If you are working with a batch template, the only available function is Save.

Follow these procedures:

- [To specify startup or shutdown methods](#)
- [To automatically update the timed SRM information](#)
- [To specify a calibration batch](#)
- [To specify device states](#)
- [To save a batch for later acquisition or processing](#)
- [To start acquisition or processing](#)
- [To view the output files](#)

❖ To specify startup or shutdown methods

1. Select a method from the System Startup Method list.

The application runs this method before running the batch. No autosampler injection takes place. This feature is not available for all instruments.

2. Select a method from the System Shutdown Method list.

The application runs this method after running the batch. This feature is not available for all instruments.

System startup method:

AS Method 1 ▼

System shutdown method:

AS Method 3 ▼

❖ To automatically update the timed SRM information

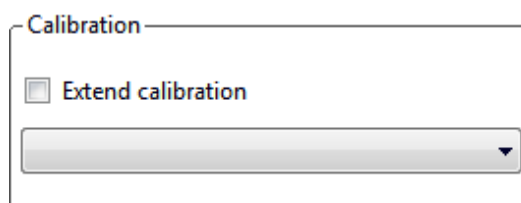
Select the **Auto TSRM Update** check box.

☐ Auto TSRM Update

When you submit the batch, the application updates the TSQ method with mass transitions, collision energy, and other appropriate data for TSRM functionality.

❖ **To specify a calibration batch**

1. In the Calibration area, select a calibration (.calx) file from the list.



Note You must acquire at least one batch with the current method to create a calibration (.calx) file.

2. To add calibration data from the current batch to the selected calibration file, select the **Extend Calibration** option.

❖ **To specify device states**

In the System Status area, select the name of the device, right-click, and then choose a device state.

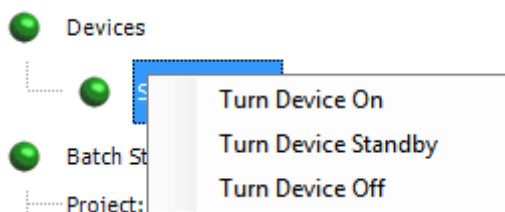






Table 10. Instrument states (Sheet 1 of 2)

Instrument state	Description
Turn Device On	Keeps the system in the On state when the current run finishes, so you can begin another run without waiting. All power and flows are maintained at operational levels. Default: On
Turn Device Standby	Keeps the system in the Standby state when the current run finishes, so you can begin another run with only a short delay between runs. Some devices do not have a Standby feature. For devices with this feature, the device enters a power-saving or consumable-saving mode, and you can switch the device back on in approximately 15 minutes. Depending on the instrument, this state turns liquid flows off but maintains heaters and other subsystems in an On state so that there is no warm-up time required when you change from Standby to On.

Table 10. Instrument states (Sheet 2 of 2)

Instrument state	Description
Turn Device Off	<p>Keeps the system in the Off state when the current run finishes. The Off state indicates that all power to the instrument, which the application can control, is turned off. This includes power to all heaters and subassemblies, but in some cases not all subassemblies.</p> <p>Some devices do not have an Off feature. For devices that do have this feature, the device enters a power-saving or consumable-saving mode, and you can switch the device back on. When several runs are queued, the application uses the system power scheme of the last submitted run.</p>
Instrument status indicators	
	Green indicates that the device is turned on or is running.
	Yellow indicates that the device is in standby mode or is waiting for contact closure.
	Red indicates that the device is turned off or that there is an error with the device.

❖ **To save a batch for later acquisition or processing**

From the Finish page, click **Save**,  **Save**.

The application saves your batch as a prepared file.

❖ **To start acquisition or processing**

1. Click **Submit**,  **Submit**.

The **Submit Options** dialog box opens.

2. To acquire (or reacquire) the submitted samples, select the **Acquire Data** check box.
 - When all submitted samples have been previously acquired, this option is (by default) not selected.
 - When one or more submitted samples have not been acquired, this option is (by default) selected.

Tip You can process the data with or without performing peak detection. For example, you might want to turn off peak detection when reprocessing samples.

3. To process the submitted samples, select the **Process Data** check box.

The application displays options for each type of method that the batch uses: Quantitation, Target Screening, Unknown Screening, or a combination of methods.

4. For each method type, select the check box for the options that you want to use.

Peak Detect: Performs peak detection for all method types. You can process the data with or without performing peak detection. For example, you might want to turn off peak detection when reprocessing samples.

Quantitate: Performs quantitation for quantitation methods.

Identify: Performs identification for unknown screening methods.

Identify and Confirm: Performs both identification and confirmation for target screening methods.

With RT Alignment: Performs retention time alignment for unknown screening methods. This produces the heat map and group averages data in the Unknown Screening View. When you select this option, the application automatically selects the Peak Detect option.

5. (Optional) Select the **Create Reports** check box.
6. (Optional with multiplexing activated) Select the **Priority Sequence** check box.

The application acquires the priority batch on the next available channel or the assigned channel.

7. (Optional without multiplexing activated) Select the **Priority Sequence** check box and then select one of the following priority options to place the batch in the queue:
 - **Next Available Batch** places the batch immediately after the currently acquiring batch.
 - **Next Available Sample** places the batch immediately after the currently acquiring sample.

Note When you select Full Sequence Submission in the Configuration console, these options are unavailable because the current batch and the current sample are, in effect, the same thing.

8. To specify the following optional parameters, click **Show Details**.
 - a. Select the **Use** check box for the device that you want to use for this acquisition.
 - b. Select the **Start Device** check box to indicate the device that will initiate communication with the other instruments.

This is usually the autosampler.
 - c. Select the **Start When Ready** check box, which starts all instruments together when they are all ready.

When you clear this check box, individual instruments can start at different times and then must wait for the last instrument to be ready.
 - d. Select the system state after it acquires the last batch: **On**, **Standby**, or **Off**.

9. To start the selected processes, click **OK**.

The selected processes begin, and the application shows the real-time display at the bottom of the current window. You can begin another batch in the Acquisition mode while you watch the real-time display of the currently acquiring batch.

IMPORTANT When your batch uses unknown screening features, after the first processing, you might see an error message stating that the total number of results is too large for a single batch (more than 500 000). Return to the Processing pages for the unknown screening features (refer to Chapter 6, “Using the Method Development Mode for Unknown Screening Methods,” in the *TraceFinder Lab Director User Guide*) and make one or more of these parameter adjustments:

- Limit the RT range.
- Shorten the signal range.
- Specify a lower value for the Number of Top Matches.
- Specify a Simple Search instead of an Exhaustive Search.
- Specify Top Peaks instead of All Peaks.
- Limit the number of search types.

Figure 13. Submit Options dialog box

Table 11. Submit Options dialog box parameters (Sheet 1 of 3)

Parameter	Description
User Name	Name of the current user.
Samples	Number of samples to be submitted for acquisition, upload, processing, or reporting.
Method Types	Lists the types of methods (quantitation, target screening, or unknown screening) used in the submitted batches.
Workflow Steps to Perform	
Acquire Data	<p>Submits the current batch to acquisition.</p> <ul style="list-style-type: none"> When all submitted samples have been previously acquired, this option is (by default) not selected. When one or more samples in the batch have not been acquired, this option is (by default) selected.

Table 11. Submit Options dialog box parameters (Sheet 2 of 3)

Parameter	Description
Process Data	<p>Processes the data for the current batch using any of the following options:</p> <p>Peak Detect: Performs peak detection for all method types. You can process the data with or without performing peak detection. For example, you might want to turn off peak detection when reprocessing samples.</p> <p>Quantitate: Performs quantitation for quantitation methods.</p> <p>Identify: Performs identification for unknown screening methods.</p> <p>Identify and Confirm: Performs both identification and confirmation for target screening methods.</p> <p>With RT Alignment: Performs retention time alignment for unknown screening methods. This produces the heat map and group averages data in the Unknown Screening View. When you select this option, the application automatically selects the Peak Detect option.</p>
Create Reports	Creates reports for the current batch.
Submit as a Priority Sequence?	
Priority Sequence	<p>With multiplexing activated, places the batch immediately after the currently acquiring batch.</p> <p>Without multiplexing activated, specifies one of the following priority options to place the batch in the queue:</p> <p>Next Available Batch: Places the batch immediately after the currently acquiring batch.</p> <p>Next Available Sample: Places the batch immediately after the currently acquiring sample.</p>
<p>Note When you select Full Sequence Submission in the Configuration console, these options are unavailable because the current batch and the current sample are, in effect, the same thing.</p>	
Acquisition pane	
Device Name	<p>Lists all configured instruments.</p> <p>If the instrument that you want to use is not configured, close the TraceFinder application, configure the instrument, and then reopen the application. You cannot configure an instrument while the TraceFinder application is running.</p> <p>Available only when you select the Acquire Data check box.</p>

Table 11. Submit Options dialog box parameters (Sheet 3 of 3)

Parameter	Description
Use	Specifies the instruments used for this acquisition. Available only when you select the Acquire Data check box.
Start Device	Specifies the instrument that initiates the communication with the other instruments. This is usually the autosampler. Available only when you select the Acquire Data check box.
Start When Ready	Starts the specified device when all the instruments are ready to acquire data. When this is cleared, individual instruments can start at different times and then must wait for the last instrument to be ready.
Post-run System State	Specifies the system state after it acquires the last batch: On (default), Standby, or Off.
Buttons	
Hide/Show Details	Collapses or expands the acquisition details of the Submit Options dialog box.
OK	Begins the selected processes.
Cancel	Closes the Submit Options dialog box without submitting any tasks.

❖ **To view the output files**

Locate the files to view from the following directories:

The application writes saved batches to the project folder:

...\\TraceFinderData\\4.0\\Projects\\...

For each acquired sample, the application writes an RSX file to the batch Data folder:

...\\TraceFinderData\\4.0\\Projects\\...\\Data

The application saves method information to the batch Methods folder:

...\\TraceFinderData\\4.0\\Projects\\...\\Methods

The application writes the reports to the batch Reports folder:

...\\TraceFinderData\\4.0\\Projects\\...*batch*\\Reports

Finish

Use the features on the Finish page to save the batch to be acquired later or acquire and process data and optionally create reports.

Figure 14. Finish page

Table 12. Finish page parameters

Parameter	Description
System Status	The System Status pane displays the following: <ul style="list-style-type: none"> Devices used for the acquisition Project, subproject, and name of the batch Number of acquired and unacquired samples in the batch Number of reports to be printed and saved as PDF, CSV, or Excel files Local method and instrument method used for the batch Number of compounds in the method
System Startup Method	The instrument methods that run before and after the batch. No autosampler injection takes place. These features are not available for all instruments.
System Shutdown Method	
Auto TSRM Update	Updates the TSQ method with mass transitions, collision energy, and other appropriate data for TSRM functionality.
Calibration	<ul style="list-style-type: none"> Use calibration: Uses the selected calibration file to process the current data. Extend calibration: Adds calibration data from the current batch to the selected calibration file.
Save	Saves the current batch as a prepared batch.
Submit	Opens the Submit Options dialog box.

Real Time Status Pane

You can access the Real Time Status pane from any mode in the application.

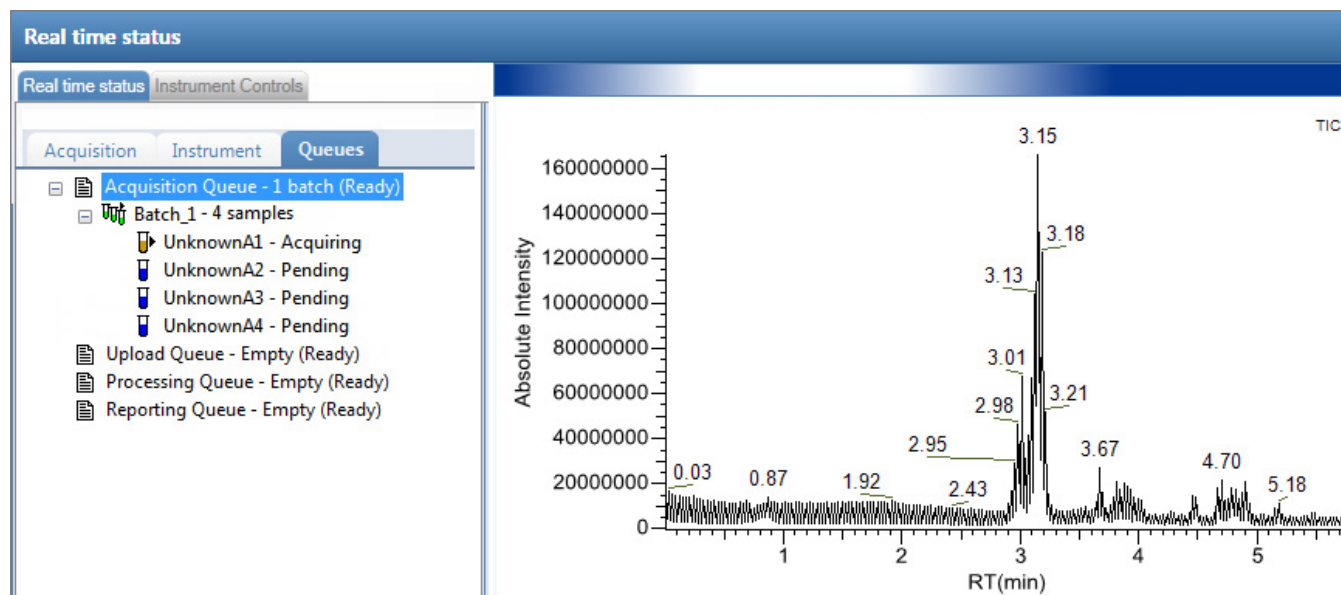
❖ **To access the Real Time Status Pane from any mode**

Click **Real Time Status** in the upper right corner of the TraceFinder window.

Real time status

The Real Time Status pane opens at the bottom of the current view.

Figure 15. Real Time Status pane



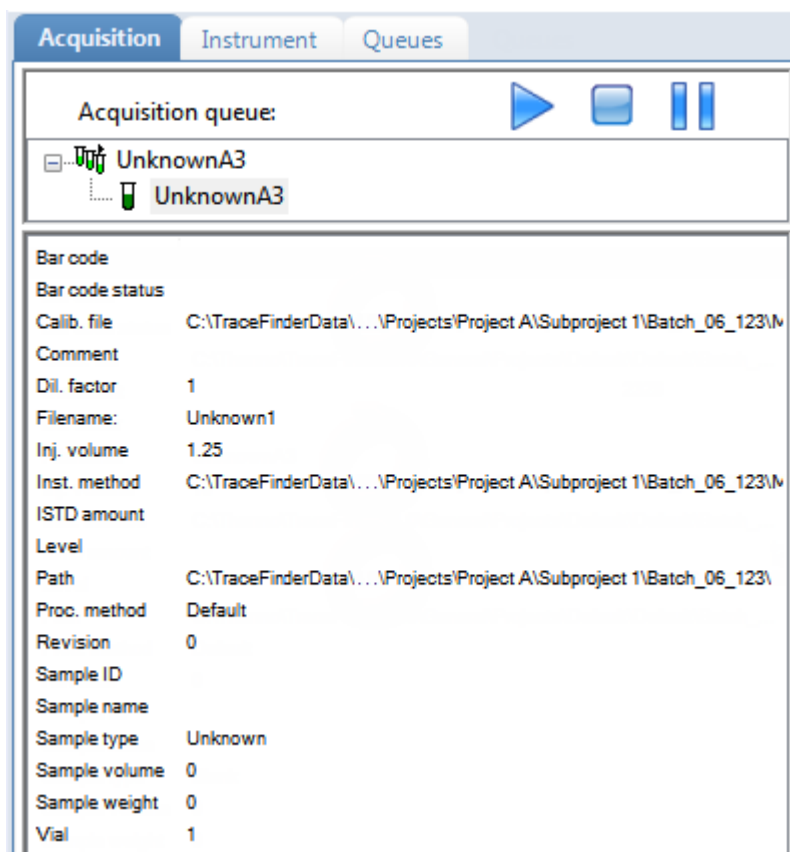
The Real Time Status pane has four pages of information and a real-time trace pane:

- [Real Time Status – Acquisition Page](#)
- [Real Time Status – Instrument Page](#)
- [Real Time Status – Queues Page](#)
- [Instrument Controls Page](#)
- [Real-Time Trace Display](#)

Real Time Status – Acquisition Page

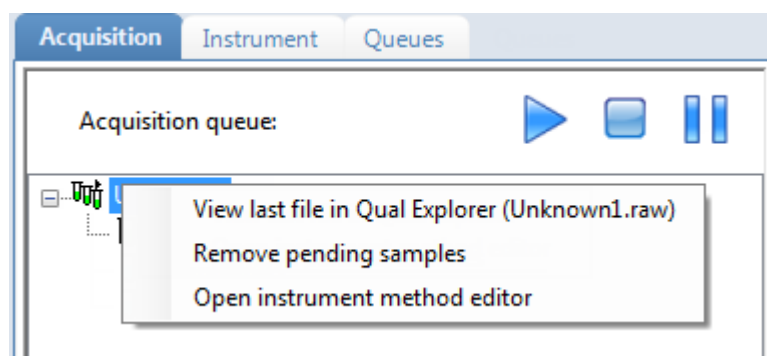
Use the Acquisition page to monitor the progress as the application acquires the samples.

Use the Start, , Stop, , or Pause, , icons to control batches in the Acquisition queue.



❖ To display the last acquired raw data file in a qualitative browser

On the Acquisition page of the real-time status pane, right-click and choose **View Last File in Qual Explorer**.



The last acquired file opens in either the FreeStyle or Qual Browser application.

❖ **To open the Instrument Setup window**

Right-click anywhere in the Acquisition page and choose **Open Instrument Method Editor**.

The Thermo Instrument Setup window opens, displaying the currently running instrument method.

For detailed information about editing instrument methods, refer to Chapter 3, “Using Instrument Methods in the Method Development Mode,” in the *TraceFinder Lab Director User Guide*.

Note Changes you make and save to the instrument method do not affect the currently running batch.

Real Time Status – Instrument Page

Use the Instrument page to monitor the currently acquiring sample.

Device Name	Status
Accela AS	Ready for Run

When you run single sample submission, this displays the sample name instead of the batch name.

❖ **To view the last acquired file in a qualitative browser**

Right-click anywhere in the top pane of the Instrument page and choose **View Last File in Qual Explorer**.

The last acquired file opens in either the FreeStyle or Qual Browser application.

❖ **To open the Instrument Setup window**

Right-click anywhere in the top pane of the Instrument page and choose **Open Instrument Method Editor**.

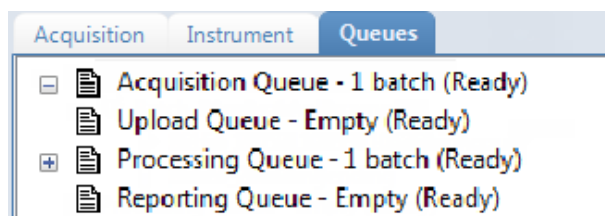
The Thermo Instrument Setup window opens, displaying the currently running instrument method.

For detailed information about editing instrument methods, refer to Chapter 3, “Using Instrument Methods in the Method Development Mode,” in the *TraceFinder Lab Director User Guide*.

Note Changes you make and save to the instrument method do not affect the currently running batch.

Real Time Status – Queues Page

Use the Queues page to monitor and control the Acquisition, Upload, Processing, and Reporting queues.



- [Queue-Level Commands](#): Pause or remove batches in any of the queues.
- [Batch-Level Commands](#): Pause or remove entire batches or samples within batches from any of the queues.
- [Additional Commands](#): Open the FreeStyle application or the Instrument Setup window.

Queue-Level Commands

Use the queue-level commands to pause or remove batches in any of the queues on the Queues page. See [Queue-Level Shortcut Menu](#).

Follow these procedures:

- [To pause all batches in a queue](#)
- [To remove a single batch from a queue](#)
- [To remove all batches in a queue](#)
- [To remove all pending batches](#)

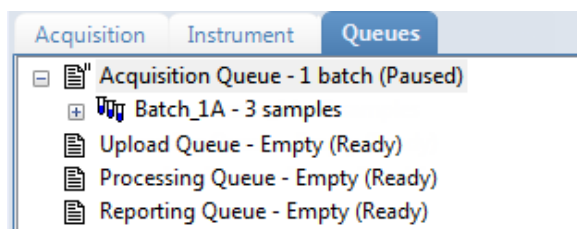
❖ **To pause all batches in a queue**

1. Select a queue (Acquisition, Upload, Processing, or Reporting).

Note When multiplexing is activated, you can have as many as four samples acquiring at once. Pausing the Acquisition queue does not affect any acquiring samples.

2. Right-click and choose **Pause Queue**.

After the current sample is complete, the application pauses all batches and samples in the specified queue. Only the selected queue is affected.



3. To restart a paused queue, select the queue, right-click, and choose **Resume Queue**.

❖ **To remove a single batch from a queue**

1. Select a queue (Acquisition, Upload, Processing, or Reporting).
2. Right-click and choose **Stop Active Batch**.

Note This command is available only when there are active batches in the queue. Paused batches and batches that contain only pending samples are not “active.”

The application confirms that you want to remove the active batch from the selected queue. After the current sample is complete, the application removes the batch and all pending samples from the queue. Only the selected queue is affected.

❖ **To remove all batches in a queue**

1. Select a queue (Acquisition, Upload, Processing, or Reporting).
2. Right-click and choose **Stop All Batches**.

The application removes all batches with pending samples from the selected queue. The current sample continues to acquire. Only the selected queue is affected.

❖ **To remove all pending batches**

1. Select a queue (Acquisition, Upload, Processing, or Reporting).
2. Right-click and choose **Remove Pending Batches**.

Note A pending batch is a batch in which all samples are pending. If any sample in the batch is active, the batch is not affected by this command.

The application removes all batches that contain only pending samples. Only the selected queue is affected.

Queue-Level Shortcut Menu

Use the commands in the shortcut menu to pause or remove batches in any of the queues on the Queues page.

Figure 16. Queue-level shortcut menu

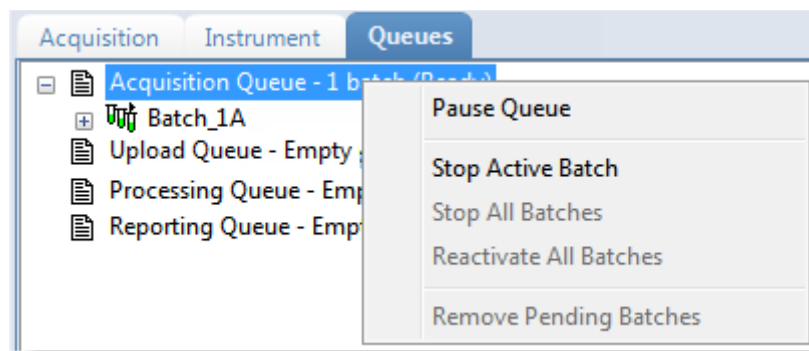


Table 13. Queue-level shortcut menu commands

Command	Description
Pause Queue	After the current sample is complete, the application pauses the specified queue. Only the selected queue is affected.
Resume Queue	Returns the paused queue to active status.
Stop Active Batch	Removes all pending samples from the specified queue. The active sample is not affected.
Stop All Batches	Removes all pending samples and batches from the specified queue. The active sample is not affected.
Reactivate All Batches	Returns all paused batches to active status.
Remove Pending Batches	Removes all pending batches from the specified queue. The active batch is not affected.

Batch-Level Commands

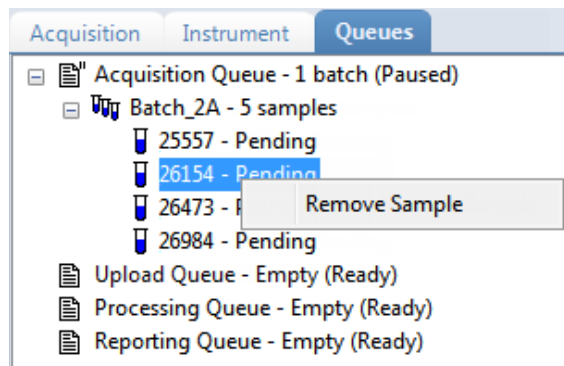
Use the batch-level commands to pause or remove entire batches or samples within batches from any of the queues on the Queues page. See [Batch-level shortcut menu](#).

Follow these procedures:

- [To remove a single pending sample from a batch](#)
- [To remove all pending samples from a batch](#)
- [To stop a batch](#)
- [To remove a pending batch](#)

❖ **To remove a single pending sample from a batch**

1. Select a pending sample.
2. Right-click the sample and choose **Remove Sample**.



The application confirms that you want to remove the selected sample from the batch and then removes the sample.

❖ **To remove all pending samples from a batch**

1. Select a batch in any of the queues (Acquisition, Upload, Processing, or Reporting).

The batch must have at least one pending sample.

2. Right-click and choose **Remove Pending Samples**.

The application confirms that you want to remove all pending samples from the batch and then removes the samples. If the batch includes only pending samples, the application removes the batch from the queue.

❖ **To stop a batch**

1. Select an active batch in any of the queues (Acquisition, Upload, Processing, or Reporting).

Note The batch must have at least one active sample and one pending sample.

2. Right-click and choose **Stop Batch**.

The application confirms that you want to remove the selected batch from the queue. After the current sample is complete, the application removes the batch and all pending samples from the queue.

❖ **To remove a pending batch**

1. Select a pending batch in any of the queues (Acquisition, Upload, Processing, or Reporting).

Note A pending batch is a batch in which all samples are pending. If any sample in the batch is active, this command is not available.

2. Right-click and choose **Remove Pending Batch**.

The application confirms that you want to remove the selected batch from the queue and then removes the batch from the queue.

Figure 17. Batch-level shortcut menu

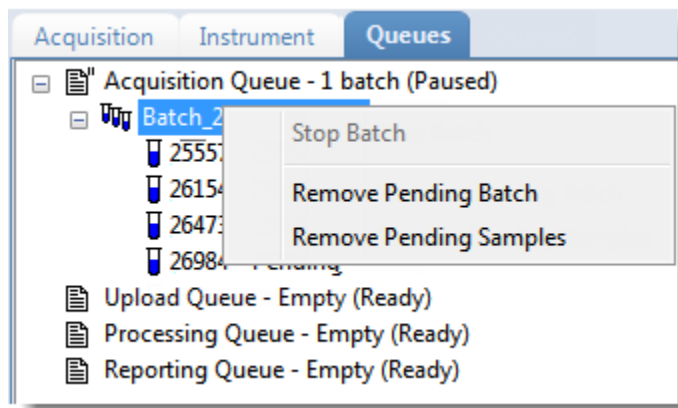


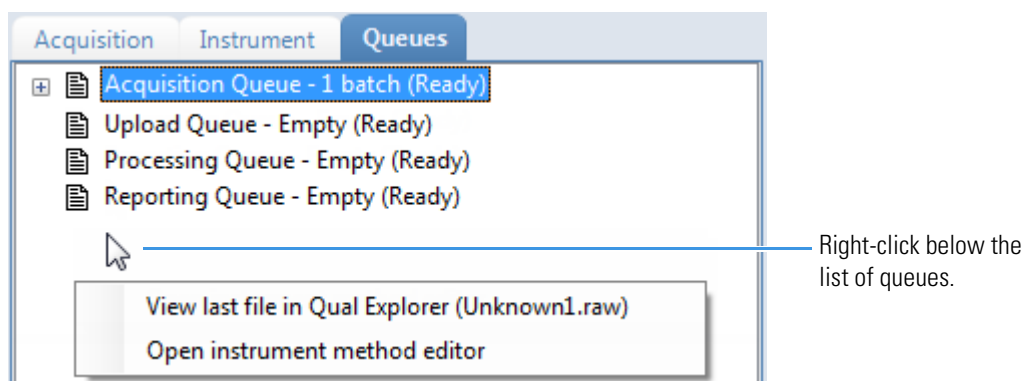
Table 14. Batch-level shortcut menu commands

Command	Description
Stop Batch	After the current sample is complete, the application removes all samples in the selected batch.
Remove Pending Batch	Removes all samples from the selected pending batch.
Remove Pending Samples	Removes all pending samples from the selected batch.

Additional Commands

You can open the FreeStyle application or the Instrument Setup window using additional commands that you access from the Queues page.

To access the shortcut menu with these commands, you must click in the white space below the list of queues. If you click on the name of or to the right of the queues, the application displays queue-, batch-, or sample-level shortcut menus.



❖ **To view the last acquired file in a qualitative browser**

Right-click below the queues list on the Queues page and choose **View Last File in Qual Explorer**.

The last acquired file opens in either the FreeStyle or Qual Browser application.

❖ **To open the Instrument Setup window**

Right-click below the queues list on the Queues page and choose **Open Instrument Method Editor**.

The Thermo Instrument Setup window opens, displaying the currently running instrument method.

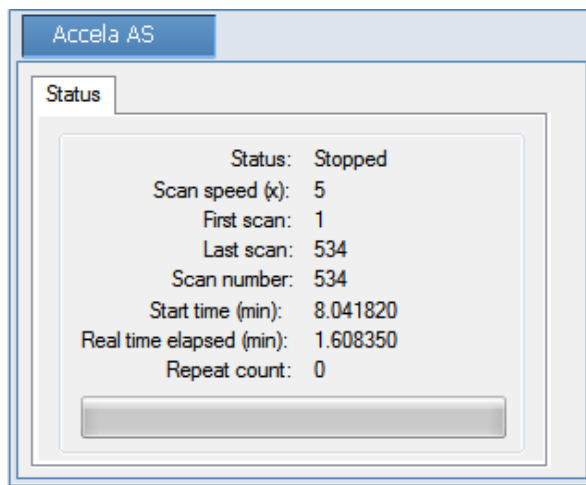
For detailed information about editing instrument methods, refer to Chapter 3, “Using Instrument Methods in the Method Development Mode,” in the *TraceFinder Lab Director User Guide*.

Note Changes you make and save to the instrument method do not affect the currently running batch.

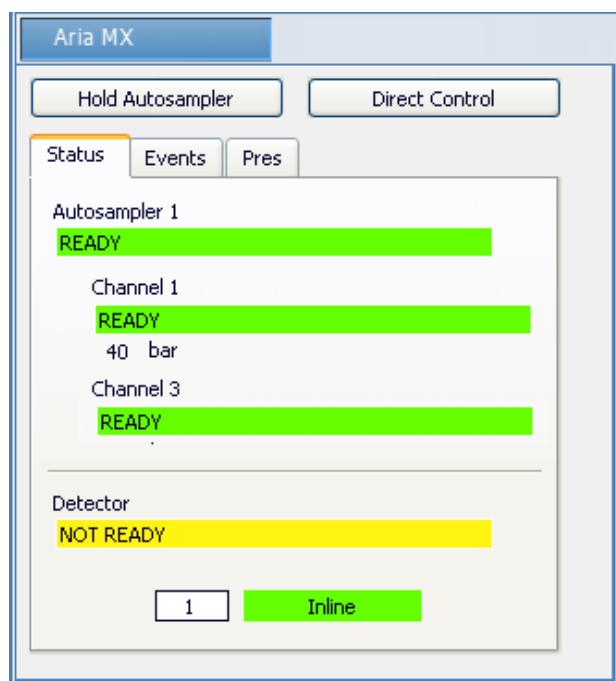
Instrument Controls Page

Use the Instrument Controls page to monitor the status of the instrument. The feedback you see on the Instrument Controls page depends on the instrument you are using. The following examples show an Accela™ autosampler and an Aria™ multiplexing device.

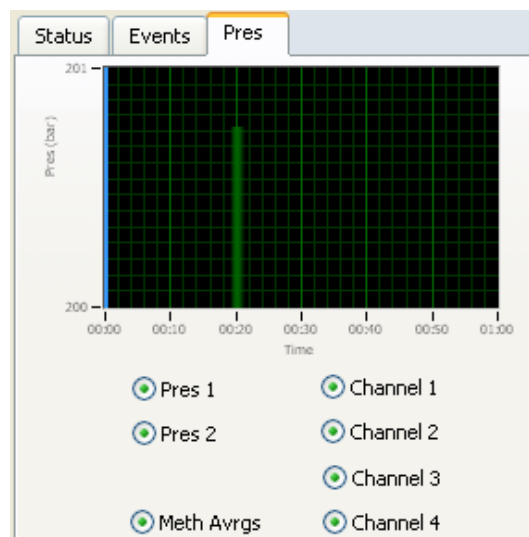
Accela Autosampler Feedback



Aria Multiplexing Feedback



Time	Ch	Msg
11:40:38.48	1	Chan Status NOT READY
11:40:36.48	1	(Channel1\Pump1) Std Pump 1
11:35:38.76		Detector NOT READY
11:35:36.30	1	Chan Status READY
11:33:24.39		System Init



Follow these procedures:

- To pause the autosampler
- To control the channels
- To open the Instrument Setup window

- To view the pressure trace
- To access the Aria multiplexing controls
- To view the last acquired file in a qualitative browser

❖ **To pause the autosampler**

1. Click **Hold Autosampler**.

The autosampler finishes the current autosampler step and then pauses. The autosampler and LC pumps continue to run.

2. To restart the autosampler, click **Hold Autosampler** again.

❖ **To control the channels**

Right-click the channel name and choose a command from the shortcut menu.

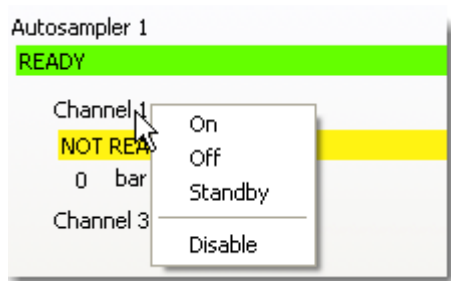


Table 15. Autosampler shortcut menu commands

Command	Description
On	Turns on a stopped pump and continues acquiring the sample list assigned to that channel.
Off	After the current sample is complete, the application stops acquiring and the pump shuts down.
Standby	After the current sample is complete, the application stops acquiring. The pump continues to run.
Disable/Enable	<p>Disable: Prevents the channel from receiving samples. When you choose Disable during a run, the application finishes the current sample on the channel and then stops.</p> <p>Enable: Allows the channel to receive samples.</p> <p>When you disable a channel that is set to On, the channel is highlighted in green and the status is READY. You can turn the channel to Off or Standby.</p>

❖ **To open the Instrument Setup window**

Right-click in the header of the Instrument Controls page and choose **Open Instrument Method Editor**.

The Thermo Instrument Setup window opens, displaying the currently running instrument method.

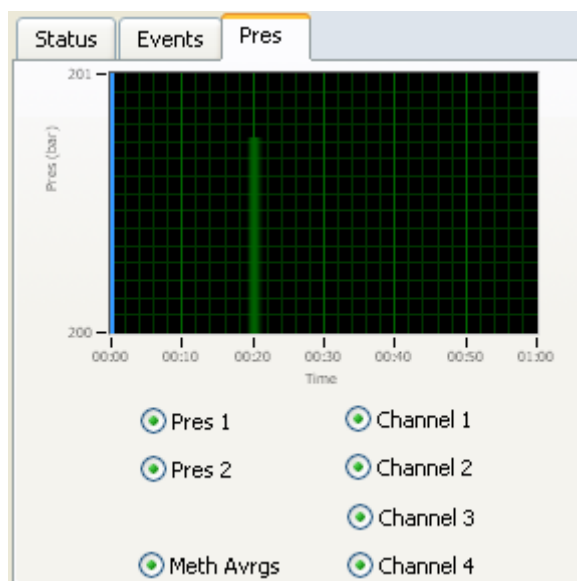
For detailed information about editing instrument methods, refer to Chapter 3, “Using Instrument Methods in the Method Development Mode,” in the *TraceFinder Lab Director User Guide*.

Note Changes you make and save to the instrument method do not affect the currently running batch.

❖ **To view the pressure trace**

1. Click the **Pres** tab.

The Pressure page displays a pump pressure graph for each sample in the batch. A fluctuation or change in the pump pressure could indicate a change in the chromatography conditions.

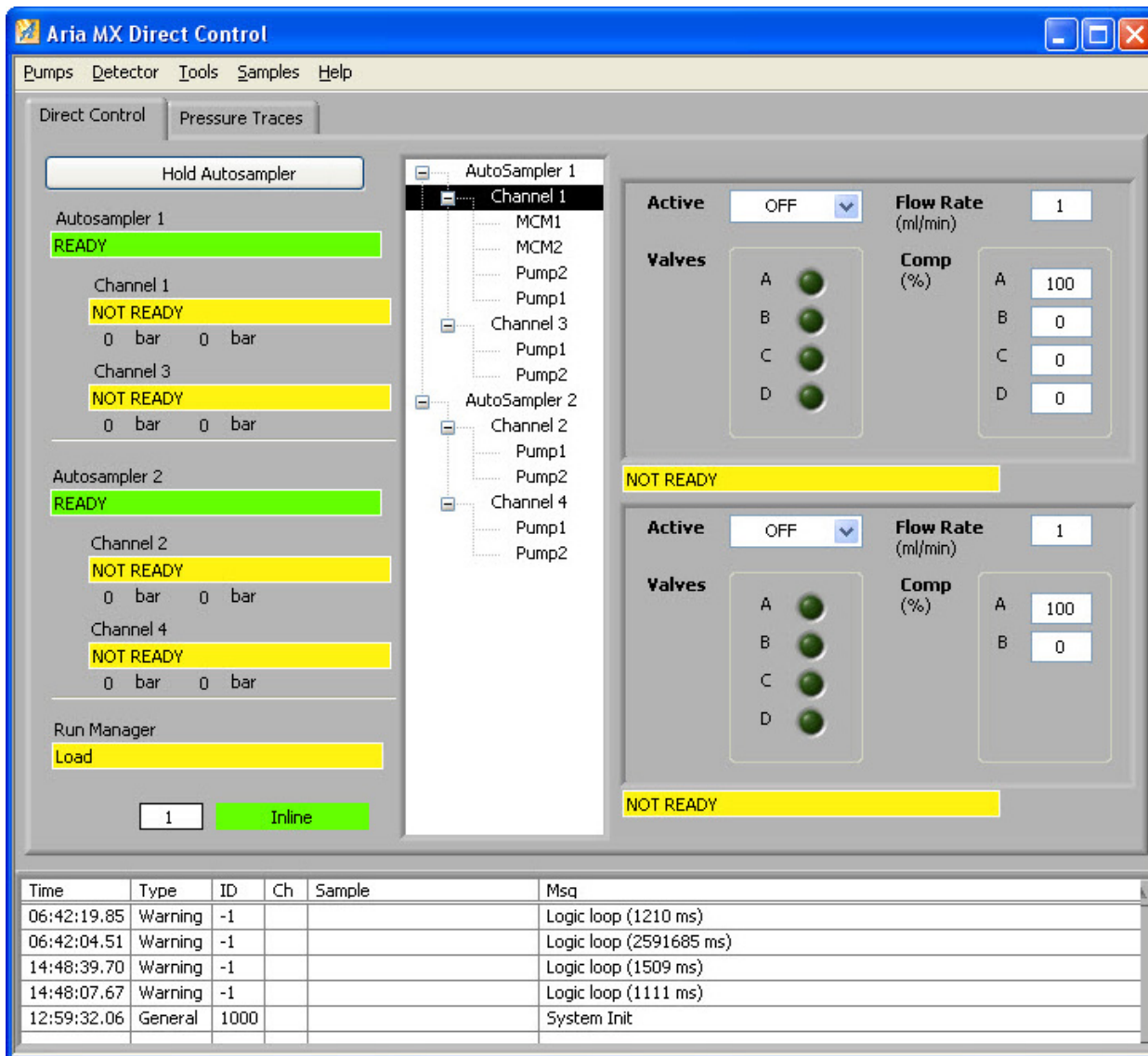


2. To view the pressure for a specific pump, select the **Pres 1** or **Pres 2** option.
By default, the pressures for all pumps are displayed.
3. To view the pressure for a specific channel, select the corresponding channel number.
By default, the pressures for all channels are displayed.

❖ **To access the Aria multiplexing controls**

Click **Direct Control**.

The Aria MX Direct Control window opens.



For detailed descriptions of the features in this window, refer to the *Transcend Systems with Xcalibur Software User Guide*.

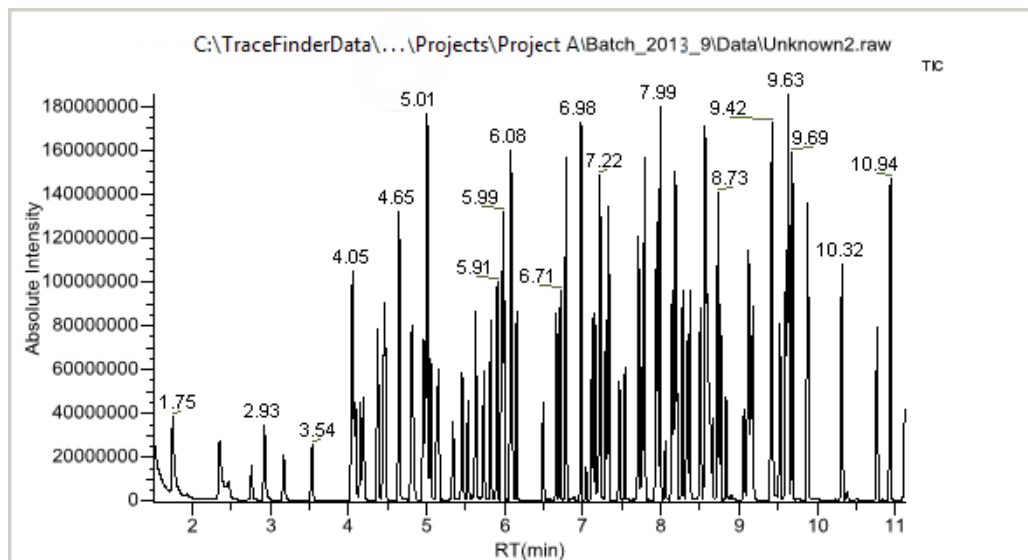
❖ **To view the last acquired file in a qualitative browser**

Right-click in the header of the Instrument Controls page and choose **View Last File in Qual Explorer**.

The last acquired file opens in either the FreeStyle or Qual Browser application.

Real-Time Trace Display

As each sample is acquired, the real-time chromatogram pane shows the retention time and intensity of the TIC trace.

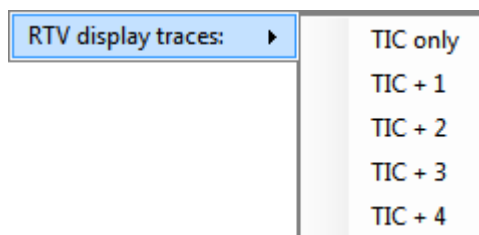


By default, the Real Time Status pane shows only the TIC trace as each sample is acquired. To observe specific traces, such as the internal standard, use the RTV Display Traces function to display multiple traces.

When you create your method, you can specify additional traces to display in the real-time viewer and in which order the traces are displayed. The application always displays the TIC trace in the top pane. See [Real Time Status Pane](#).

❖ To display multiple traces

Right-click the chromatogram pane and choose the number of traces to display.

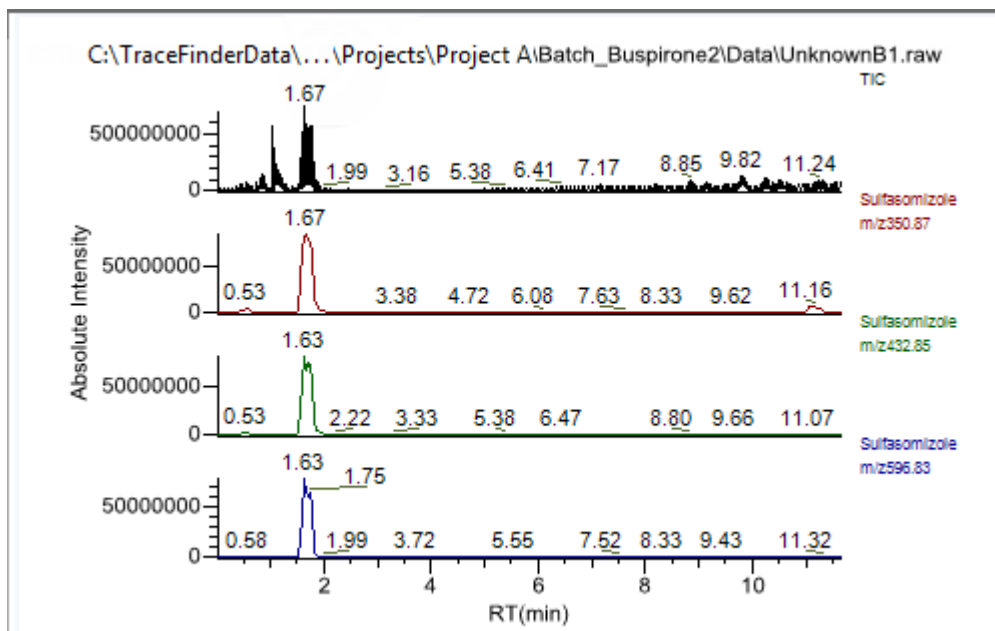


The chromatogram pane displays real-time chromatograms for the selected number of traces.

The TIC is always displayed at the top. When there are more traces than can fit in the pane, you can scroll through the traces.

For each trace, the application displays the mass or precursor mass. See [Real-time trace display with multiple traces](#).

Figure 18. Real-time trace display with multiple traces



Sample Types

The TraceFinder application uses the following sample types in all sample definitions and reports.

Table 16. Sample type definitions

Sample type	Definition
Negative	Contains no target compounds but might contain an ISTD when you use the internal standard quantitative analysis technique. By analyzing a blank sample, you can confirm that there are no residual compounds in the solvent system that can cause erroneous results.
Unextracted	Similar to a Negative sample, but contains target compounds. By analyzing a sample, you can confirm that there are no residual compounds in the solvent system that can cause erroneous results.
Calibrator	(Calibration standard) Contains known amounts of all target compounds. The purpose of a standard is to measure the response of the instrument to the target compounds so that the processing application can generate a calibration curve for each compound.
QC	(Quality Check) Contains a known amount of one or more specific target compounds. The application places check standard samples in the sequence so that it can test quantitative analysis results for quality assurance purposes. After the application analyzes the QC sample, it compares the measured quantity with the expected value and an acceptability range. The quantitative analysis of a QC sample is classified as <i>passed</i> if the difference between the observed and expected quantities is within the user-defined tolerance. A QC sample is classified as <i>failed</i> if the difference between the observed and expected quantities is outside the user-defined tolerance.
Solvent	Contains only solvent.
Specimen	Used for the quantitative analysis of samples.
Hydrolysis	Checks the degradation of compounds dissolved in water.

3 Using the Acquisition Mode

Sample Types

Using the Analysis Mode for Quantitation Batches

Use the features of the Analysis mode to do the following:

- Create quantitation batches.
- Submit quantitation batches for acquisition, processing, or report generation.
- Review quantitation batches, batch data, reports, and local methods.

IMPORTANT TraceFinder 4.1 uses the same data as TraceFinder 4.0. By default, the application stores the method, project, template, and compound database data for the 4.1 release in the TraceFinderData\4.0 folder.

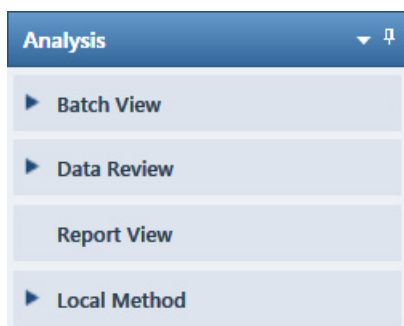
Contents

- [Working in the Batch View for Quantitation Batches](#)
- [Working in Data Review for Quantitation Batches](#)
- [Working in the Report View for Quantitation Batches](#)
- [Working in the Local Method View for Quantitation Batches](#)

❖ To access the Analysis mode

Click **Analysis** in the navigation pane.

The Analysis navigation pane opens.



Working in the Batch View for Quantitation Batches

In the Batch View, you can manually create and edit a new quantitation batch or open and edit a previously saved batch. When you submit a batch, you can acquire and process data and optionally create reports for the submitted samples.

The Analysis mode includes a toolbar:



Use the [Toolbar](#) or the equivalent commands in the [Batch View Shortcut Menu](#) to create the sample list and submit samples for acquisition.

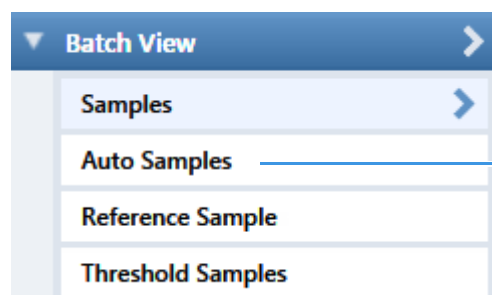
See the following topics:

- [Samples Page for Quantitation Batches](#)
- [Auto Samples Page for Quantitation Batches](#)
- [Reference Sample Page for Quantitation Batches](#)
- [Threshold Samples Page for Quantitation Batches](#)

❖ To open the Batch View

1. Click **Analysis** in the navigation pane of the current mode.
2. Click **Batch View**.

The Batch View navigation pane opens.



Available only when you activate Intelligent Sequencing in the Configuration console

Samples Page for Quantitation Batches

To open the Samples page, click **Samples** in the Batch View navigation pane.

For details about the Samples page, see the following topics:

- [Samples Page Features](#)
- [Creating a New Quantitation Batch](#)
- [Editing a Quantitation Batch](#)
- [Submitting a Quantitation Batch](#)

Samples Page Features

The Samples page is divided into two panes:

- Samples pane
Use the [Samples Pane](#) to create a batch.
- Compound Active Status pane
Use the [Compound Active Status Pane](#) to make specific compounds active or inactive.

Samples pane

The screenshot shows the 'Batch View - Batch_A [Quan]' window. At the top, there is a 'Local Method' dropdown set to 'Method_Quan', an 'Update' button, an 'Instrument' field set to 'Thermo Scientific Instrument', and a 'User' field. Below this is a table with 15 columns: Status, Filename, Sample type, Groups, Qual Processing, Level, Sample ID, Sample name, Vial position, Injection volume, Conversion Factor, Channel, Barcode Expected, Barcode Actual, and Sample Volume. The table contains three rows of data for 'Unknown1', 'Unknown2', and 'Unknown3'. Below the table is a 'Compound Active Status' section with a sub-table containing two rows: one for 'FENTHION-CE20-R20-TL75-QED' at RT 1.91, and another for 'Sulfisomidine' at RT 2.72. Both are marked as 'Active' with checked checkboxes.

Compound Active Status pane

Tip To resize the panes, drag the separator that divides the panes.

Samples Pane

The samples pane includes the following features:

- [Column Display](#)
- [Status Indicators](#)
- [Groups](#)
- [Qual Processing](#)
- [Sample Weight Calculation](#)
- [Instrument Methods](#)
- [Toolbar](#)
- [Batch View Sample List](#)
- [Batch View Shortcut Menu](#)

Column Display

The sample list contains many columns of information. You can scroll to see all the columns, and you can customize which ones to display and their display order.

❖ To scroll the sample list

Use the horizontal scroll bar at the bottom of the sample list to view all the information.

When you use the scroll bar at the bottom of the sample list, the following columns remain fixed: Status, Filename, Sample Type, Groups, Qual Processing, Level, Sample ID, and Sample Name, while the other columns scroll right and left.

❖ To customize the column display

1. Right-click the sample list and choose **Modify Columns**.

The **Modify Columns dialog box** opens.

2. Use the arrow buttons to move all the columns that you want displayed to the Displayed Columns pane.

These columns appear to the right of the sample list's fixed columns: Status, Filename, Sample Type, Groups, Qual Processing, Level, Sample ID, and Sample Name.

3. To arrange the order of the columns, do the following:

- a. In the Displayed Columns pane, select a column name.
- b. Use **Up** or **Down** to move the selected column up or down in the list.

The first column in the list represents the leftmost column in the Batch View sample list, and the last column in the list represents the rightmost column in the Batch View sample list.

Note The following columns are fixed: Status, Filename, Sample Type, Groups, Qual Processing, Level, Sample ID, and Sample Name.

4. To change the width of a column, do the following:

- a. In the Displayed Columns pane, select the column width.

5	Sample ID	100
▶ 6	Sample name	100
7	Vial position	100

- b. Type a new value for the width.

5. Repeat [step 4](#) for all columns whose widths you want to change, and click **OK**.

The columns in the sample list immediately reflect your changes. The application uses these settings for all sample lists in the Batch View.

Figure 19. Modify Columns dialog box

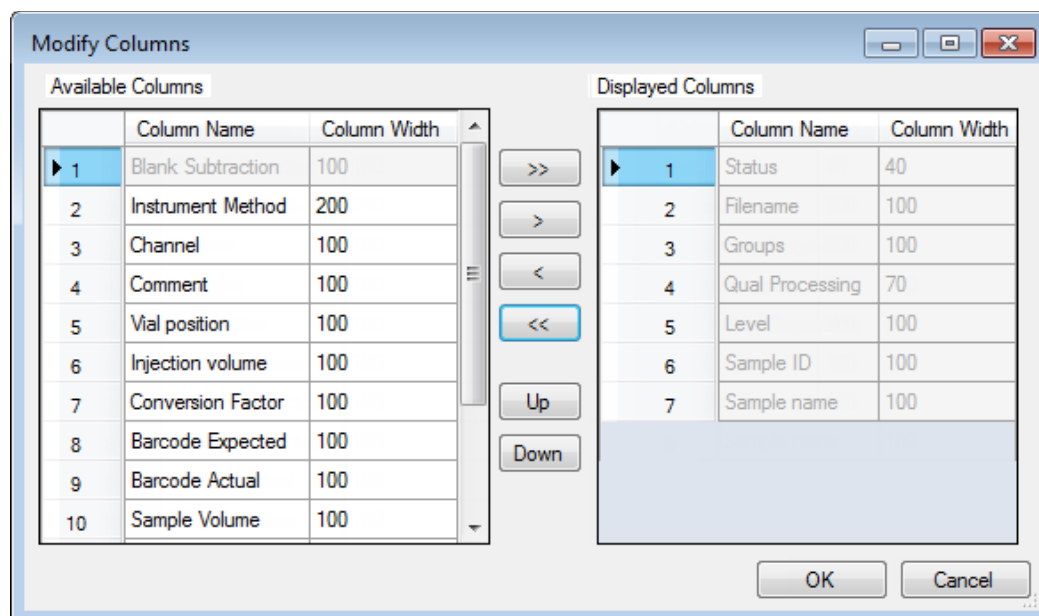




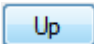
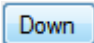






Table 17. Button descriptions for the Modify Columns dialog box




Button	Description
	Moves all columns to the Displayed Columns pane.
	Moves the selected column to the Displayed Columns pane.
The following buttons apply to all columns, except for those that are fixed: Status, Filename, Sample Type, Groups, Qual Processing, Level, Sample ID, and Sample Name.	
	Moves the selected column to the Available Columns pane.
	Moves all columns except fixed columns.
	Moves the selected column name in the Displayed Columns pane one row up in the column order.
	Moves the selected column name in the Displayed Columns pane one row down in the column order.

Status Indicators

Status indicators show the current status of each sample during the acquisition and processing.

-  Sample is not acquired.
-  Sample is acquired but not processed.
-  Sample is acquired and processed.
-  Sample is currently acquiring.

Local Method: Method_Alprazolam Update

	Status	Filename	Sample type	Groups	Qual Processing	Sample ID	Sample name	V p
1		Solvent_A	Solvent		<input type="checkbox"/>			1
2		Cal_5	Calibrator		<input type="checkbox"/>			2
3		Unknown1	Specimen		<input type="checkbox"/>			3

Status indicators




Groups

Use the Groups feature to assign samples to a group.

❖ To create a group



- For each sample, type the name of a group in the Groups column.
Repeat this for each sample that you want to include in a group.
- Create as many groups as you want.

Note To assign a sample to multiple groups, separate the groups with a comma.

Status	Filename	Groups
	Benzo26473	groupB, groupA
	Benzo25557	groupB
	Benzo26154	groupB, groupA

Qual Processing

Use the Qual Processing feature to indicate samples to be processed with the qualitative peak processing criteria specified in the method.

Status	Filename	Sample type	Groups	Qual Processing
	Benzo25557	Specimen		<input checked="" type="checkbox"/>
	Benzo25558	Specimen		<input checked="" type="checkbox"/>

The Qualitative View displays processed data for the selected samples. See [Qualitative View for Quantitation Batches](#).

Sample Weight Calculation

Use the sample weight features to calculate the conversion factor for a sample. The application uses different methods to calculate the conversion factor for liquid or solid calculation types.

Liquid: $SampleVolume \div DilutionFactor$

Solid: $(SampleVolume \times DilutionFactor) \div SampleWeight$

Manual: The application does not calculate the Conversion Factor. Instead, you can enter the Conversion Factor value.

Follow these procedures:

- [To display the features for calculating sample weight](#)
- [To calculate the conversion factor for a liquid sample](#)
- [To calculate the conversion factor for a solid sample](#)
- [To manually specify the conversion factor for a sample](#)

❖ To display the features for calculating sample weight

If the Conversion Factor, Sample Volume, Dilution Factor, Sample Weight, Calculation Type, and Final Units columns are not visible, right-click and choose **Enable Sample Weight Calculation**.

Conversion Factor	Sample Volume	Dilution Factor	Sample Weight	Calculation Type	Final Units
1.000	1	1	1	Liquid	
1.000	1	1	1	Solid	
1.000	1	1	1	Manual	

❖ To calculate the conversion factor for a liquid sample

Note The application uses the following formula to calculate the Conversion Factor:
 $SampleVolume \div DilutionFactor$

1. From the **Calculation Type** list, select **Liquid**.

For a liquid sample, the Sample Weight value is not editable.

2. In the Sample Volume column, type the volume in ng/mL for your sample.
3. In the Dilution Factor column, type the value for the dilution.

For example, if you have 1000 ng/mL of a substance that is too concentrated for the mass spectrometer, you can dilute it by 1000. Then your injection volume is 1, your conversion factor is 1000, and your sample amount is 1000.

4. In the Final Units column, type the units that you want to use for the calculated amount in the Data Review view or in reports.

❖ To calculate the conversion factor for a solid sample

Note The application uses the following formula to calculate the Conversion Factor:
 $(SampleVolume \times DilutionFactor) \div SampleWeight$

1. From the **Calculation Type** list, select **Solid**.
2. In the Sample Weight column, type the weight in ng for your sample.
3. In the Sample Volume column, type the volume in ng/ml for your sample.
4. In the Dilution Factor column, type the value for the dilution.

For example, if you have 1000 ng/ml of a substance that is too concentrated for the mass spectrometer, you can dilute it by 1000. Then your injection volume is 1, your conversion factor is 1000, and your sample amount is 1000.

5. In the Final Units column, type the units that you want to use for the calculated amount in the Data Review view or in reports.

❖ To manually specify the conversion factor for a sample

Note The application uses the specified conversion factor when it calculates the amount for the sample.

1. From the **Calculation Type** list, select **Manual**.

For a manually calculated sample, the only available columns are the Conversion Factor and the Final Units.


2. In the Conversion Factor column, type the conversion factor to use for your sample.
3. In the Final Units column, type the units that you want to use for the calculated amount in the Data Review view or in reports.

Instrument Methods

Use the Instrument Methods column to specify instrument methods for the samples.

Note By default, the Instrument Method column is not displayed in the Batch View sample list.




❖ To specify instrument methods for samples

1. Display the Instrument Method column in the sample list:
 - a. Right-click the sample list and choose **Modify Columns**.
The Modify Columns dialog box opens.
 - b. In the Available Columns pane, select **Instrument Method**.
 - c. Click  to move the Instrument Method column to the Displayed Columns pane.
 - d. Click **OK**.

The application displays the Instrument Method column, defaulting to the instrument method specified in the master method.

2. Click the Instrument Method column and select an instrument method from the list.
This list contains all the available instrument methods. Instrument methods from external sources are prefixed with “Ext:”.

You can specify a different instrument method for each sample.

Batch View - Batch_8266 [Quan]						
Local Method: Method_Quan		Update		Instrument: Thermo Scientific Instrument		
	Status	Groups	Filename	Sample type	Qual Processing	Instrument Method
4			Unknown4	Specimen	<input type="checkbox"/>	Ext: Instrument1
5			Unknown5	Specimen	<input type="checkbox"/>	Ext: Instrument2
6			Unknown6	Specimen	<input type="checkbox"/>	Ext: Instrument3

When you submit the batch for acquisition, the application saves a copy of the selected instrument methods to the following folders:

External instrument methods:

...\\TraceFinderData\\4.0\\Projects\\...\\batch\\Methods\\method\\ExternalMethods

Local instrument methods:

...\\TraceFinderData\\4.0\\Projects\\...\\batch\\Methods\\method

Toolbar

The Analysis mode includes this toolbar for creating and submitting a batch.



Table 18. Toolbar icons

Icon	Description
	Adds the specified number of new, empty samples to the end of the sample list. See the instructions To add samples to the list .
	Inserts a new, empty sample or samples above the selected sample. See the instructions To insert samples into the list .
	Removes the selected samples from the sample list. See the instructions To remove samples from the list .
	Adds imported samples from a CSV, an XML, or an SLD file to the sample list. See the instructions To import samples into the list .
	Submits only the selected samples for acquisition, processing, or report generation. See the instructions To submit samples in the batch .
	Submits the batch for acquisition, processing, or report generation. See the instructions To submit samples in the batch .
	Submits only the selected peaks for processing. See the instructions To submit selected peaks for processing .
	Opens the Acquisition mode where you can use a batch template to define a standard sequence composed of various sample types to be assembled into a batch of samples. See Working in Data Review for Quantitation Batches .
	Opens the Acquisition mode where you can create a batch template that contains the basic settings and sample types for your batches. See Using the Acquisition Mode .
	Opens the Quick Acquisition window where you can quickly submit a single sample. See Appendix A, "Using Quick Acquisition."
	Opens the Audit Viewer where you can view audit logs. See Chapter 8, "Using the Audit Viewer." Available only when you enable Auditing in the Administrator Console. Refer to the instructions in the <i>TraceFinder Administrator Console User Guide</i> .

Batch View Sample List

The sample list displays all the quantitative data for the samples of a batch.

Status indicators for each sample indicate if the sample is currently acquiring, not acquired, acquired, or processed.

The sample list includes the following columns of information:

Figure 20. Batch View sample list

<div> Local Method: Method_Benzos_Quan Update Instrument: Thermo Scientific Instrument </div>										
	Status	Filename	Sample type	Groups	Qual Processing	Level	Sample ID	Sample name	Vial position	Injection volume
1		UnknownA1	Specimen		<input type="checkbox"/>		1		CStk1-01:7	10.0
2		UnknownA2	Specimen		<input type="checkbox"/>		1		CStk1-01:7	10.0
3		UnknownA3	Specimen		<input type="checkbox"/>		1		CStk1-01:7	10.0

Calculation Type	Conversion Factor	Dilution Factor	Sample Weight	Sample Volume	Final Units
Liquid ▼	1.000	1	1	1	
Liquid ▼	1.000	1	1	1	
Liquid ▼	1.000	1	1	1	

Instrument Method	Channel	Barcode Expected	Barcode Actual	Comment
Instrument1 ▼	Auto			
Instrument1 ▼	Auto			
Instrument1 ▼	Auto			

Table 19. Batch View sample list columns (Sheet 1 of 2)





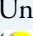
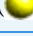

Column	Description
Status	<p> Sample is not acquired.</p> <p> Sample is acquired but not processed.</p> <p> Sample is acquired and processed.</p> <p> Sample is currently acquiring.</p> <p>Note When you include unknown screening features in the quantitation method and you choose to process with only the quantitation criteria, the Sample View shows the Status for the samples as acquired and processed () , and the Unknown Screening View shows the Status for the samples as acquired but not processed () .</p> <p>Note When you include unknown screening features in the quantitation method and you choose to process with only the unknown screening criteria, both the Sample View and the Unknown Screening View show the Status for the samples as acquired and processed () .</p>
Groups	Threshold group to which a sample belongs. Samples can be viewed by group in the Comparative View of Data Review.
Filename	Name of the raw data file that contains the sample data.
Sample Type	<p>Defines how the application processes the sample data. Each sample is classified as one of the following sample types: Specimen, QC, Solvent, Calibrator, Hydrolysis, Unextracted, or Negative.</p> <p>Default: Specimen</p>
Qual Processing	Indicates samples to be processed with the qualitative peak processing criteria specified in the method. The Qualitative View displays processed data for the selected samples.
Level	The level defined for a calibration sample or quality control sample.
Sample ID	A user-defined, alphanumeric string that identifies a sample.
Sample Name	A user-defined name that identifies a sample.
Vial Position	The tray vial number used for an autosampler acquisition.
Injection Volume	<p>The injection volume (in microliters) of the injected sample.</p> <p>When you are using an autosampler, you can set the default injection volume in the Autosampler dialog box in the Instrument View. The minimum and maximum injection volumes that you can use depend on the autosampler you configure. The usable range depends on the injection mode and might be smaller than the displayed range.</p> <p>The Injection Volume value set in the master method overwrites the value in the instrument method.</p> <p>Valid range: 0.1 through 5000 µL</p>

Table 19. Batch View sample list columns (Sheet 2 of 2)

Column	Description
Calculation Type	<p>Liquid: The application calculates the Conversion Factor as</p> $\text{SampleVolume} \div \text{DilutionFactor}$ <p>Solid: The application calculates the Conversion Factor as</p> $(\text{SampleVolume} \times \text{DilutionFactor}) \div \text{SampleWeight}$ <p>Manual: Sample Volume, Dilution Factor, Sample Weight, and Final Units columns are not available, and the Conversion Factor value is editable.</p>
Conversion Factor	<p>Editable only when Calculation Type is Manual.</p> <p>Default: 1</p>
Sample Volume	Default: 1
Dilution Factor	Default: 1
Sample Weight	<p>Available only when Calculation Type is Solid.</p> <p>Default: 1</p>
Final Units	<p>Specifies the calculated amount in the Data Review view or in reports.</p> <p>Default: 1</p>
Instrument Method	Specifies the instrument to use for the acquisition. This column is hidden by default. To display this column, see To customize the column display .
Channel	Specifies the channel on which the sample was run. If the sample is not acquired, the value is Pending. The Channel column is available only when you have activated multiplexing in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the <i>TraceFinder Lab Director User Guide</i> .
Barcode Expected	A user-entered barcode for the vial.
Barcode Actual	An actual barcode for the vial. This value is not editable.
Comment	A user-defined comment for the sample.


Batch View Shortcut Menu

The Batch View includes a shortcut menu for creating a batch.

Table 20. Batch View shortcut menu commands (Sheet 1 of 2)

Command	Description
Add Sample	Adds a single empty row to the sample grid.
Insert Sample	Inserts a single empty row to the sample grid above the selected row.
Insert Copy Sample	Copies the currently selected row and inserts a copy above the row.
Reinject Selected Samples	Creates a copy of the selected sample and appends INJ001 to the file name. Additional reinjections of the same sample are numbered INJ002, INJ003, and so forth.
Remove Selected Samples	Removes selected samples from the sample grid.
Import Samples	Opens the Sample Import Tool. See To import samples into the list .
Browse in Raw File (Move)	Opens a dialog box where you can select a raw data file to use for the selected sample row. The application removes the raw data file from the source location.
Browse in Raw File (Copy)	Opens a dialog box where you can select a raw data file to use for the selected sample row. The application copies the raw data file from the source location.
Map Raw Files to Samples	Opens a dialog box where you can select multiple raw data files to use for the selected sample rows.
Copy Down	Copies the value in the selected row to all rows below it. This command is available only when you have selected a value that can be copied down.
Fill Down	Enters sequential values in the column starting with the value in the selected row and ending with the last row in the column. This command is available only when you have selected a value that can be filled down.
Modify Columns	Opens the Modify Columns Dialog Box .
Hide Sample Weight Calculation	Displays or hides the Sample Volume, Dilution Factor, Sample Weight, Calculation Type, and Final Units columns.
Copy	Copies the data in the selected rows or columns to the Clipboard. Use this command to copy sample information into a text editor or spreadsheet application. You cannot paste this data back into the Batch View sample list.

Table 20. Batch View shortcut menu commands (Sheet 2 of 2)

Command	Description
Copy with Headers	<p>Copies the data in the selected rows or columns and the associated column headers to the Clipboard. Use this command to copy sample information into another text editor or spreadsheet application. You cannot paste this data back into the sample list.</p> <p>For example</p>  <p>Copy With Headers from TraceFinder</p> <p>Paste into Excel spreadsheet</p>
Paste	Pastes a single column of copied data from another text editor or spreadsheet application into the selected column.
Undo Last Paste	Removes the last pasted item in the Batch View.
Export to CSV File	Opens the Save As dialog box where you can save the current sample list to a CSV file.
Edit Instrument Method	<p>Opens the Instrument Setup window where you can edit the parameters of the instrument method.</p> <ul style="list-style-type: none"> When you edit an external method, the application updates the method in the ...\\Xcalibur\\methods folder. When you edit an internal method, the application updates the method in the ...\\TraceFinderData\\4.0\\Projects\\...\\batch\\Methods\\method folder. <p>For detailed information about editing instrument methods, refer to Chapter 3, “Using Instrument Methods in the Method Development Mode,” in the <i>TraceFinder Lab Director User Guide</i>.</p>
View Sample in Qual Explorer	Displays the selected sample in the qualitative explorer application that you configured as your default in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the <i>TraceFinder Lab Director User Guide</i> .

Compound Active Status Pane

In the Compound Active Status pane, you can choose specific compounds to be active or inactive.

❖ To set a compound as active or inactive

1. In the sample list, select a sample.

All compounds in the selected sample are listed in the Compound Active Status pane.

Compound Active Status

	RT	Compound	Active
▶ 1	1.91	FENTHION-CE20-R20-TL75-QED	<input checked="" type="checkbox"/>
2	2.72	Sulfisomidine	<input checked="" type="checkbox"/>

The default active/inactive status is determined by the identification settings in the local method. For information about setting the identification parameters, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

- To display compounds alphabetically, right-click and choose **Sort by Compound Name**.
 - To display compounds from shorter to longer retention time, right-click and choose **Sort by Retention Time**.
2. Select or clear the **Active** check box for the compound.

To change the active/inactive status in the Data Review view, see [Inactive and Excluded Compounds](#).

Compound Active Status



You can specify which compounds are active or inactive in the Local Method View or the Batch View.

Figure 21. Active and inactive compounds in the Local Method View

Local Method View - Batch_1_Method_Benzos					
Master method: Method_Benzos					
Acquisition List	Identification	Detection	Calibration	Calibration levels	QC Check levels
	RT	Compound	Compound type	Active	CAS No
1	1.91	FENTHION-CE20-R20-TL...	Target Compound	<input type="checkbox"/>	55389
▶ 2	2.72	Sulfisomidine	Target Compound	<input checked="" type="checkbox"/>	515640

For details about setting the status on the Identification page, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

Figure 22. Active and inactive compounds in the Batch View

Batch View - Batch_2 [Quan]								
Local Method: Method_Benzo1_Quan				Update				
	Status	Filename	Sample type	Groups	Qual Processing	Level	Sample ID	ne
▶ 1		Benzo25555	Specimen		<input type="checkbox"/>		CStk1-01:3	
2		Benzo25557	Specimen		<input type="checkbox"/>		CStk1-01:4	

Compound Active Status			
	RT	Compound	Active
▶ 1	5.88	1-Butanol, 4-(butylnitrosoami...	<input type="checkbox"/>
2	16.68	3,3'-Diaminodiphenylmethane	<input checked="" type="checkbox"/>
3	2.91	4-Methyl-2,6,7-trioxa-1-phosp...	<input checked="" type="checkbox"/>

For details about setting the status in the Batch View, see [Compound Active Status Pane](#).

Creating a New Quantitation Batch

In the Batch View, you can create a new batch.

Follow these procedures:

- [To create a new batch](#)
- [To add samples to the list](#)
- [To insert samples into the list](#)
- [To import samples into the list](#)
- [To remove samples from the list](#)
- [To copy a sample](#)
- [To reinject a sample](#)
- [To edit sample values](#)
- [To browse in raw data files](#)
- [To customize the column display](#)

❖ To create a new batch

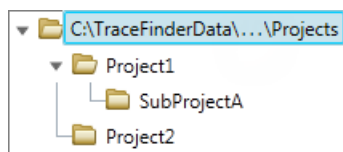
1. Choose **File > New > Batch** from the main menu.

The [Create New Batch Dialog Box](#) opens, displaying all drives that contain projects.

4 Using the Analysis Mode for Quantitation Batches

Working in the Batch View for Quantitation Batches

2. Select a drive from the list.



Tip The application displays all configured and enabled repositories.

3. Select the folder where you want to store your batch.

Tip To activate the Create button, you must enter a unique batch name. If the Create button is not activated, you have entered a batch name that is already used.

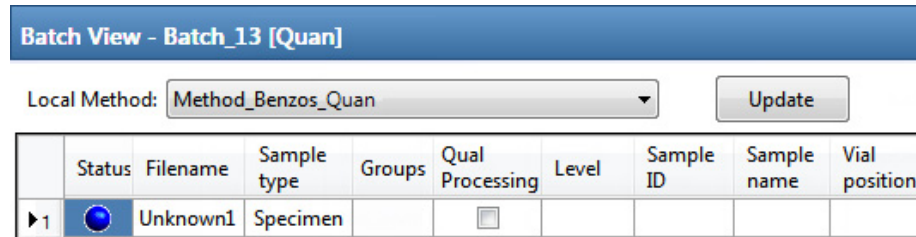
To create a new folder for the storage location, see [Editing Folders for Batches](#).

4. Select **Quan** from the Type list.

The batch list displays all batches in the selected folder. The Method list displays all methods for the quantitation type.

5. Select a master method from the Master Method list.
6. Click **Create**.

A new batch opens with one Specimen sample.




The batch name in the title bar indicates that you are creating a quantitation batch.

When you create a batch using a quantitation method that includes unknown screening features, the title bar indicates that the method includes unknown screening.




❖ To add samples to the list

1. To add a single sample row, right-click the sample list and choose **Add Sample**.
2. To add multiple sample rows, select the number of rows and then click the **Add Sample** icon, .





The application adds the specified number of new, empty samples to the end of the sample list.

❖ To insert samples into the list

Select the sample above which you will insert new, Specimen samples, and then do one of the following:


- To insert a single sample row, right-click and choose **Insert Sample**.
- To insert multiple sample rows, select the number of rows and then click the **Insert Sample** icon .

The application inserts the Specimen samples above the selected sample.

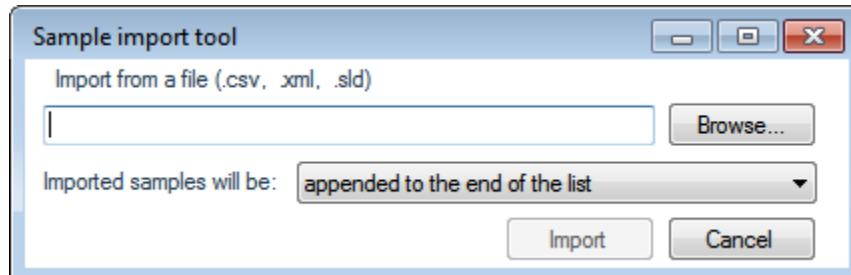
	Status	Filename	Sample type	Groups
1		cal_std_5	Calibrator	
2		Unknown2	Specimen	
3		Unknown1	Specimen	
4		cal_std_10	Calibrator	

Inserted samples

❖ To import samples into the list

1. Choose **Batch > Import Samples** from the main menu, or click the **Import Samples** icon, .

The Sample Import Tool dialog box opens.



From this dialog box, you can import samples from a CSV, an XML, or an SLD file.

2. Click **Browse** and select a CSV, an XML, or an SLD file that contains the samples to import.
3. From the Imported Samples Will Be list, select **Appended to the End of the List** or **Inserted at the Selected Row**.
4. Click **Import**.

The Sample Import Tool dialog box closes, and the application adds the specified samples to the sample list.

4 Using the Analysis Mode for Quantitation Batches

Working in the Batch View for Quantitation Batches

When you import samples from an Xcalibur sequence file (.sld), the application makes the following column name substitutions.

Xcalibur column	TraceFinder column
Position	Vial Position
Inj Vol	Injection Volume
Dil Factor	Conversion Factor

When you import samples from an Xcalibur sequence file (.sld), the application makes the following sample type substitutions.

Xcalibur sample type	TraceFinder sample type
Blank	Negative
Std Bracket	Calibrator

5. (Optional) When using multiplexing, select a channel for each imported sample.

Imported samples default to Auto.

Note The Channel column is available only when you have activated multiplexing in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.

❖ To remove samples from the list

1. Select the samples that you want to remove.

Tip Use the CTRL or SHIFT keys to select multiple samples.

2. Right-click and choose **Remove Selected Samples**.

❖ To copy a sample

1. Select the sample that you want to copy.
2. Right-click and choose **Insert Copy Sample**.

The application inserts the copy above the selected sample.

❖ To reinject a sample

1. In the sample list, select the sample that you want to reinject.
2. Right-click and choose **Reinject This Sample**.

The application creates a copy of the selected sample and appends INJ001 to the file name. Additional reinjections of the sample are numbered INJ002, INJ003, and so forth. The application copies all parameter values from the original sample.

❖ **To edit sample values**

1. For each sample, do one of the following:

Type a new file name over the current filename.

–or–

Double-click the Filename column and locate a raw data file to use for the sample.

–or–

Right-click and choose **Browse in Raw File**, and then locate a raw data file to use for the sample.

By default, the application sets the Sample Type to Unknown.

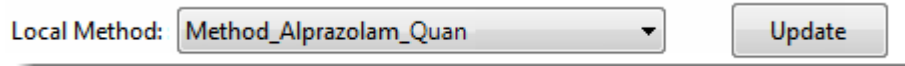
2. For each sample, click the Sample Type column and select a sample type from the list.

Available sample types			
Specimen	Hydrolysis	Solvent	QC
Unextracted	Calibrator	Negative	

3. For each Calibrator or QC sample, select a level from the Level list.

The sample levels are defined in the master method. If there are no levels to select in the Level list, do the following:

- a. Return to the Method Development mode.
- b. Open the method.
- c. Click **Compounds** in the navigation pane.
- d. Click the **Calibration Levels** tab.
- e. Add the levels.
- f. Save the method.
- g. Return to the Batch View in the Analysis mode, and then click **Update**.



The application updates the local method with the new sample levels.

For detailed instructions about specifying calibration levels, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

4. Type a vial position in the Vial Position column for each sample.
5. Type a volume in the Injection Volume column for each sample.

The minimum injection volume value allowed is 0.1 µL; the maximum injection volume value allowed is 5000 µL.

6. (Optional) Type or edit the values for the remaining columns.

Note When you use the scroll bar at the bottom of the sample list, the following columns remain fixed: Status, Filename, Sample Type, Groups, Qual Processing, Level, Sample ID, and Sample Name, while the other columns scroll right and left.

To automatically copy or fill values in these columns, see [Appendix B, “Using Copy Down and Fill Down.”](#)

❖ To browse in raw data files

1. Do one of the following:

Double-click the Filename column.


—or—

Right-click and choose **Browse in Raw File**.

The What Raw File Would You Like to Use dialog box opens.

2. Select a raw data file to use for the sample or use the CTRL key to select multiple files, and then click **Open**.

The application overwrites the selected, unacquired sample in the batch with the first “browsed in” file and adds any additional browsed in files below the selected sample.

For all browsed-in raw data files, the application sets the Status to Acquired, , and sets the Sample Type to Specimen.

Note You cannot overwrite an acquired sample. When you select a sample that is acquired, the application adds all browsed in files below the selected sample.

❖ To customize the column display

1. Right-click the Batch View sample list and choose **Modify Columns**.

The [Modify Columns dialog box](#) opens.

2. Use the arrow buttons to move all the columns that you want displayed to the Displayed Columns pane.

All the columns you select appear after the Status, Filename, Sample Type, Groups, Qual Processing, Level, Sample ID, and Sample Name columns.

3. To arrange the order of the columns, do the following:
 - a. In the Displayed Columns pane, select a column name.
 - b. Use **Up** or **Down** to move the selected column up or down in the list.

The first column in the list represents the leftmost column in the Batch View sample list, and the last column in the list represents the rightmost column in the Batch View sample list.

Note The following columns are fixed: Status, Filename, Sample Type, Groups, Qual Processing, Level, Sample ID, and Sample Name.

4. To change the width of a column, do the following:
 - a. In the Displayed Columns pane, select the column width.

7	Vial position	69
▶ 8	Injection volume	100
9	Calculation Type	100

- b. Type a new value for the width.
5. When you have completed your changes, click **OK**.

The columns in the sample list immediately reflect your changes. The application uses these settings for all sample lists in the Batch View.

Modify Columns Dialog Box

Use the Modify Columns dialog box to select the columns that you want to display.

Figure 23. Modify Columns dialog box

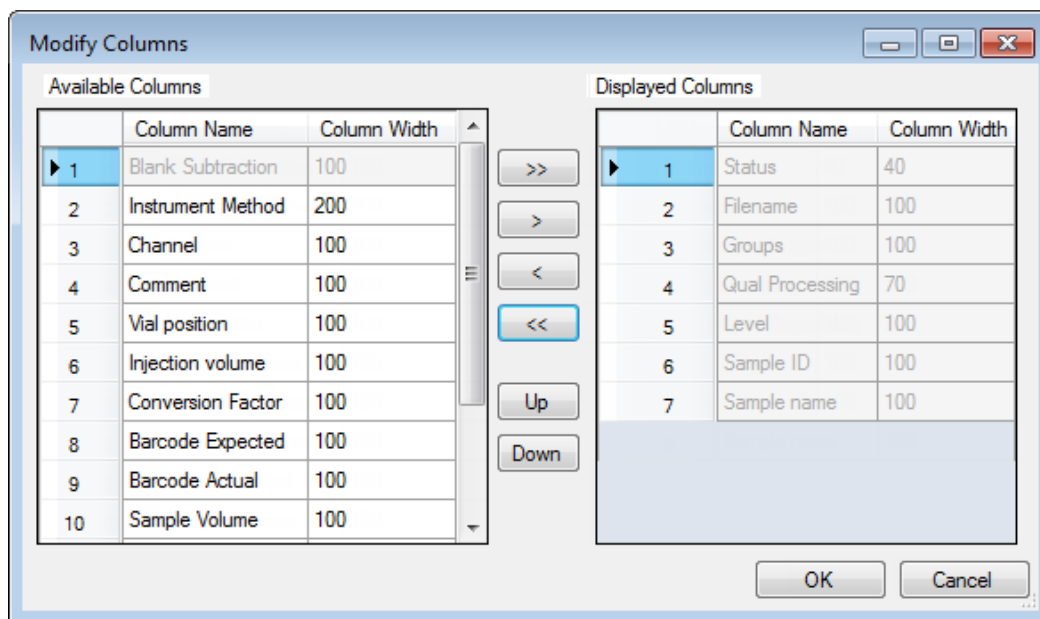




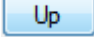
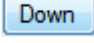
Table 21. Button descriptions for the Modify Columns dialog box (Sheet 1 of 2)

Button	Description
>>	Moves all columns to the Displayed Columns pane.
>	Moves the selected column to the Displayed Columns pane.

4 Using the Analysis Mode for Quantitation Batches

Working in the Batch View for Quantitation Batches

Table 21. Button descriptions for the Modify Columns dialog box (Sheet 2 of 2)

Button	Description
The following buttons apply to all columns, except for those that are fixed: Status, Filename, Sample Type, Groups, Qual Processing, Level, Sample ID, and Sample Name.	
	Moves the selected column to the Available Columns pane.
	Moves all columns except those that are fixed.
	Moves the selected column name in the Displayed Columns pane one row up in the column order.
	Moves the selected column name in the Displayed Columns pane one row down in the column order.

Create New Batch Dialog Box

Use the Create New Batch dialog box to select a folder and method for your batch and to name the new batch.

Figure 24. Create New Batch dialog box

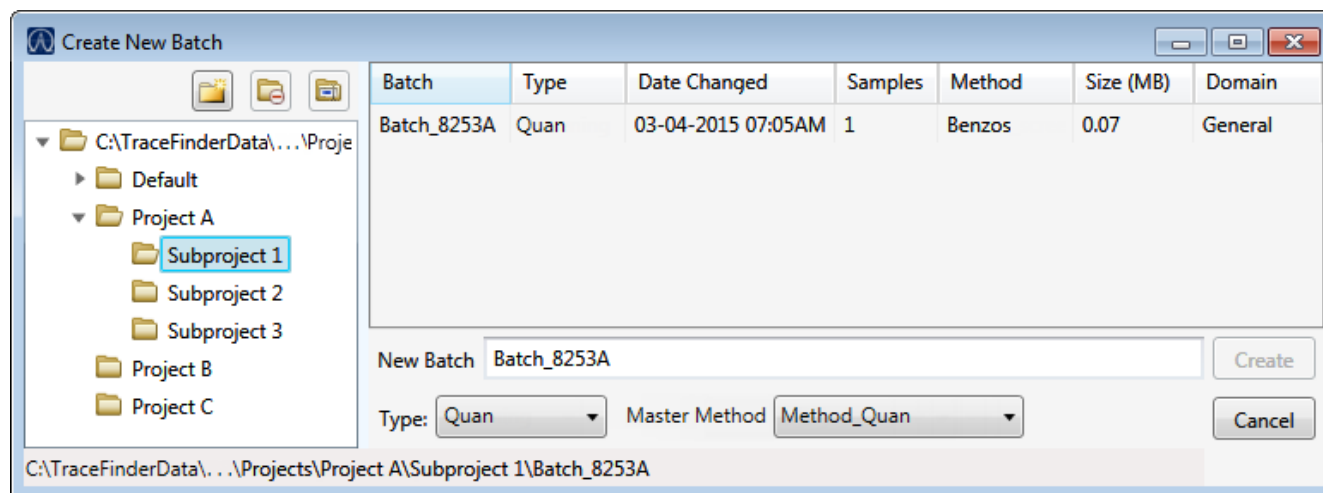


Table 22. Create New Batch dialog box parameters (Sheet 1 of 2)




Parameter	Description
 Create New Folder	<p>Adds one of the following:</p> <ul style="list-style-type: none">• When a drive is selected, adds a new project-level folder to the drive.• When a project folder is selected, adds a subproject-level folder to the selected project.• When a subproject folder is selected, adds a lower-level folder to the subproject. <p>Or, you can right-click and choose Create Folder.</p>

Table 22. Create New Batch dialog box parameters (Sheet 2 of 2)

Parameter	Description
	With no confirmation prompt, immediately removes the selected folder.
Delete Folder	You cannot delete a folder that contains lower-level folders; you must delete the lower-level folders first. Or, you can right-click and choose Delete.
	Renames the selected folder.
Rename Folder	Or, you can right-click and choose Rename.
Batch table	
Batch	Name of batches in the selected project.
Type	Type of batch: Quan, Screening, or Unknown Only.
Date Changed	Date that the batch was last updated.
Samples	Number of samples in the batch.
Method	Name of the method used to create the batch.
Size	Size of the batch in megabytes.
Domain	TraceFinder domain in which the batch was created.
New batch parameters	
New Batch	Name of the new batch to create. Note If the Create button is not activated, you have entered a name that is already used or you have not selected a method.
Type	Type of batch to create: Quan, Screening, or Unknown Only.
Method	Method used to create the new batch.
Path	Path to the project in the TraceFinderData\4.0\Projects folder where the batch is created.
Buttons	
Create	Creates the specified batch and opens the Batch View for the new batch.
Cancel	Closes the Create New Batch dialog box without creating a batch.


Editing Folders for Batches

From the Create New Batch dialog box, you can create new folders for your batches. You can also delete or rename folders.

Use these procedures:

- [To create new project folders](#)
- [To delete project folders](#)
- [To rename project folders](#)

❖ To create new project folders

1. In the Create New Batch dialog box, select the folder for which you will create a new lower-level folder.
 - You can select the main TraceFinderData\4.0\Projects folder and create a new folder under it.
 - You can select one of the existing folders and create a lower-level folder under it.
2. Click the **Create Folder** icon, .


The application adds a new lower-level folder to the selected folder.

3. Select the new folder name and type a name for the folder.

Folder names are limited to 30 characters and can contain spaces and special characters, except for the following special characters: \ / : + ? " < >

Note After you add a lower-level folder, you cannot rename the parent folder.


❖ To delete project folders

1. In the Create New Batch dialog box, select the folder to delete.
2. Click the **Delete Folder** icon, .

With no confirmation prompt, the application immediately removes the selected folder.

Note You cannot delete folders that contains lower-level folders; you must delete the lower-level folders first.

❖ To rename project folders

1. In the Create New Batch dialog box, select the folder to rename.
2. Click the **Rename Folder** icon, .

Note You cannot rename folders that contain lower-level folders.

3. Type a new name for the folder and press ENTER.

The application saves the new folder name.

Editing a Quantitation Batch

In the Batch View, you can open a saved batch and edit the sample list. You can add samples, edit samples, or remove samples. If the batch has already been acquired, you can select specific samples for reinjection. If the batch has unacquired samples when you complete your edits, you can save it as a “ready to acquire” batch.

Follow these procedures:

- To open a saved batch
- To open a recent batch
- To edit samples in a batch
- To reinject a sample from a previously acquired batch

❖ **To open a saved batch**

1. Choose **File > Open > Batch** from the main menu.

The **Open Batch dialog box** opens.

2. Select a project and a subproject.
3. Select **Quan** or **Any** from the Type list.

The batch list displays all batches created with quantitation methods (or all methods of all types when you select Any).

4. Select a batch from the list.
5. Click **Open**.

The selected batch opens in the Batch View.

Open Batch Dialog Box

Use the Open Batch dialog box to select a batch to open.

Figure 25. Open Batch dialog box

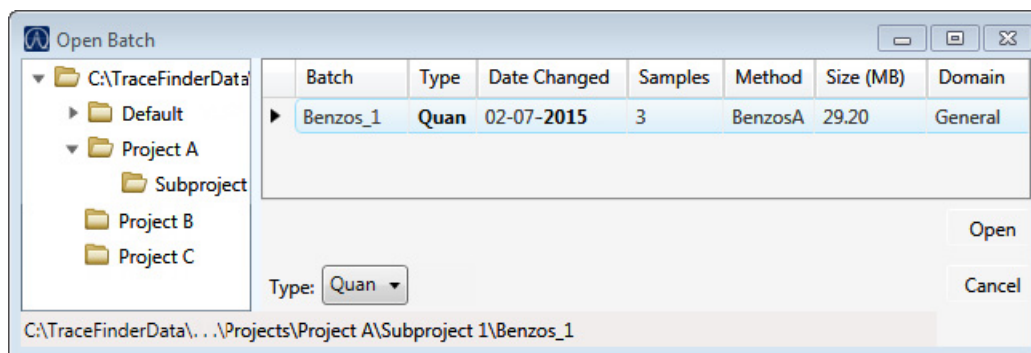


Table 23. Open Batch dialog box parameters (Sheet 1 of 2)

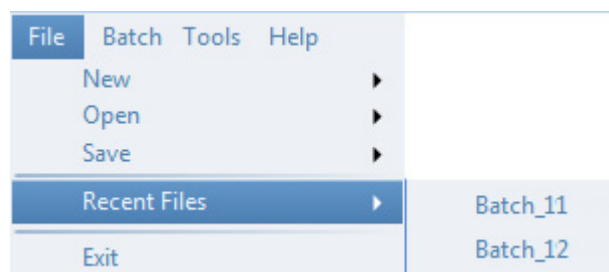
Parameter	Description
Batch	Name of batches in the selected project.
Type	Type of batch: Quan, Screening, or Unknown Only.
Date Changed	Date the batch was last updated.

Table 23. Open Batch dialog box parameters (Sheet 2 of 2)

Parameter	Description
Samples	Number of samples in the batch.
Method	Name of the method used to create the batch.
Size	Size of the batch in megabytes.
Domain	TraceFinder domain in which the batch was created.
Path	Path to the project in the TraceFinderData\4.0\Projects folder where the batch is stored.
Buttons	
Type	Type of batch to display in the Batch list: Quan, Screening, Unknown Only, or Any.
Open	Opens the Batch View for the selected batch.
Cancel	Closes the Open Batch dialog box without opening a batch.

❖ To open a recent batch

Choose **File > Recent Files > *batch*** from the main menu.



The selected batch opens in the Batch View.

❖ To edit samples in a batch

Use the commands described in [Working in the Batch View for Quantitation Batches](#).





You can add new samples, edit samples, or delete samples.

❖ To reinject a sample from a previously acquired batch

1. In the sample list, select the sample that you want to reinject.
2. Right-click and choose **Reinject This Sample**.

The application creates a copy of the selected sample and appends INJ001 to the file name. Additional reinjections of the sample are numbered INJ002, INJ003, and so forth. The application copies all parameter values from the original sample.

A green status icon indicates previously acquired samples (acquired and processed), and the sample name is grayed out. A blue status icon indicates samples created for reinjection (not acquired).

	cal_50_INJ001	Calibrator	10
	cal_50	Calibrator	10
	cal_10_INJ001	Calibrator	10
	cal_10	Calibrator	10

When you submit all samples in this batch, the application acquires all samples (including previously acquired samples).

- To save this batch with the new samples for reinjection, choose **File > Save > Batch** from the main menu.

The batch is saved as a prepared batch that is ready to submit. You can open this batch from the Reinject Samples page in the Acquisition mode and submit the batch. The application acquires only the samples that have not been previously acquired.

Submitting a Quantitation Batch


In the Batch View, you can submit an entire batch or only selected samples in the batch. When you submit a batch for acquisition and processing, you can choose to create reports for the submitted samples. See [Submit Options dialog box](#).

For a description of commands in the shortcut menu, see [Batch View shortcut menu commands](#).

Follow these procedures:



- [To submit selected peaks for processing](#)
- [To submit samples in the batch](#)
- [To view the output files](#)

❖ To submit selected peaks for processing

- In the Peak List, select the **Selected** check box for the peaks that you want to process.
- Click the **Submit ... for Processing** icon, .

The application processes the selected peaks and updates the data in the Data Review panes.

❖ To submit samples in the batch

- Do one of the following:
 - To submit all samples in the batch, click the **Submit Batch** icon, .
 - To submit specific samples, select the samples and click the **Submit Selected Samples** icon, .

The [Submit Options dialog box](#) opens.

Note You can also submit only selected peaks for processing. See [To submit selected peaks for processing](#).

2. To acquire (or reacquire) the submitted samples, select the **Acquire Data** check box.
 - When all submitted samples have been previously acquired, this option is (by default) not selected.
 - When one or more samples in the batch have not been acquired, this option is (by default) selected.

Tip You can process the data with or without performing peak detection. For example, you might want to turn off peak detection when reprocessing samples.

3. To process the submitted samples, select the **Process Data** check box.

The application displays processing options for the method.

If the quantitation method includes unknown screening features, the application also displays unknown screening options. For specifying processing data for batches that also use unknown screening features, see [To submit samples in the batch](#) of Chapter 6, “Using the Analysis Mode for Unknown Screening Batches.”

4. Select the check box for the quantitation options that you want to use.

Peak Detect: Performs peak detection. You can process the data with or without performing peak detection. For example, you might want to turn off peak detection when reprocessing samples.

Quantitate: Performs quantitation.

5. (Optional) Select the **Create Reports** check box.
6. (Optional with multiplexing activated) Select the **Priority Sequence** check box.

The application acquires the priority batch on the next available channel or the assigned channel.

7. (Optional without multiplexing activated) Select the **Priority Sequence** check box and then select one of the following priority options to place the batch in the queue:
 - **Next Available Batch** places the batch immediately after the currently acquiring batch.
 - **Next Available Sample** places the batch immediately after the currently acquiring sample.

Note When you select Full Sequence Submission in the Configuration console, these options are unavailable because the current batch and the current sample are, in effect, the same thing.

8. To specify the following optional parameters, click **Show Details**.

- a. Select the **Use** check box for the device that you want to use for this acquisition.
- b. Select the **Start Device** check box to indicate the device that will initiate communication with the other instruments.

This is usually the autosampler.

- c. Select the **Start When Ready** check box, which starts all instruments together when they are all ready.

When this is cleared, individual instruments can start at different times and then have to wait for the last instrument to be ready.

- d. Select the system state after it acquires the last batch: **On**, **Standby**, or **Off**.

9. To start the selected processes, click **OK**.

The selected processes begin, and the application shows the real-time display at the bottom of the current window. You can begin another batch in the Analysis mode while you watch the real-time display of the currently acquiring batch.

IMPORTANT When your batch also uses unknown screening features, after the first processing, you might see an error message stating that the total number of results is too large for a single batch (more than 500 000). Return to the Processing pages for the unknown screening features (refer to Chapter 6, “Using the Method Development Mode for Unknown Screening Methods,” in the *TraceFinder Lab Director User Guide*) and make one or more of these parameter adjustments:

- Limit the RT range.
- Shorten the signal range.
- Specify a lower value for the Number of Top Matches.
- Specify a Simple Search instead of an Exhaustive Search.
- Specify Top Peaks instead of All Peaks.
- Limit the number of search types.

Submit Options Dialog Box

Use the Submit Options dialog box to submit a batch for acquisition, processing, or reporting.

Figure 26. Submit Options dialog box

Table 24. Submit Options dialog box parameters (Sheet 1 of 3)

Parameter	Description
User Name	Name of the current user.
Samples	Number of samples to be submitted for acquisition, processing, or reporting.
Workflow Steps to Perform	
Acquire Data	<p>Submits the current batch to acquisition.</p> <ul style="list-style-type: none"> When all submitted samples have been previously acquired, this option is (by default) not selected. When one or more samples in the batch have not been acquired, this option is (by default) selected.

Table 24. Submit Options dialog box parameters (Sheet 2 of 3)

Parameter	Description
Process Data	<p>Processes the data for the current batch using any of the following options:</p> <p>Peak Detect: Performs peak detection. You can process the data with or without performing peak detection. For example, you might want to turn off peak detection when reprocessing samples.</p> <p>Quantitate: Performs quantitation.</p> <p>Note Quantitation methods can include unknown screening features. For specifying processing data for batches that also use unknown screening features, see To submit samples in the batch.</p>
Create Reports	Creates reports for the current batch.
Submit as a Priority Sequence?	
Priority Sequence	<p>With multiplexing activated, places the batch immediately after the currently acquiring batch.</p> <p>Without multiplexing activated, specifies one of the following priority options to place the batch in the queue:</p> <p>Next Available Batch: Places the batch immediately after the currently acquiring batch.</p> <p>Next Available Sample: Places the batch immediately after the currently acquiring sample.</p> <p>Note When you select Full Sequence Submission in the Configuration console, these options are unavailable because the current batch and the current sample are, in effect, the same thing.</p>
Acquisition pane	
Device Name	<p>Lists all configured instruments.</p> <p>If the instrument that you want to use is not configured, close the TraceFinder application, configure the instrument, and then reopen the application. You cannot configure an instrument while the TraceFinder application is running.</p> <p>Available only when you select the Acquire Data check box.</p>
Use	<p>Specifies the instruments used for this acquisition.</p> <p>Available only when you select the Acquire Data check box.</p>
Start Device	<p>Specifies the instrument that initiates the communication with the other instruments. This is usually the autosampler.</p> <p>Available only when you select the Acquire Data check box.</p>

Table 24. Submit Options dialog box parameters (Sheet 3 of 3)

Parameter	Description
Start When Ready	Starts the specified device when all the instruments are ready to acquire data. When this is cleared, individual instruments can start at different times and then must wait for the last instrument to be ready.
Post-run System State	Specifies the system state after it acquires the last batch: On (default), Standby, or Off.
Buttons	
Hide/Show Details	Collapses or expands the acquisition details of the Submit Options dialog box.
OK	Begins the selected processes.
Cancel	Closes the Submit Options dialog box without submitting any tasks.

❖ To view the output files

Locate the files to view from the following directories:

The application writes saved batches to the project folder:

...\\TraceFinderData\\4.0\\Projects\\...

For each acquired sample, the application writes an RSX file to the batch Data folder:

...\\TraceFinderData\\4.0\\Projects\\...\\Data

The application saves method information to the batch Methods folder:

...\\TraceFinderData\\4.0\\Projects\\...\\Methods

The application writes the reports to the batch Reports folder:

...\\TraceFinderData\\4.0\\Projects\\...*batch*\\Reports

Saving a Batch to a New Location

You can move the current batch to a different project folder, or you can make a copy of the current batch and save the copy to a different project folder.

Follow these procedures:

- [To save a batch to another project folder](#)
- [To move a batch to another folder](#)
- [To create a new project folder](#)
- [To delete project folders](#)
- [To rename project folders](#)

❖ To save a batch to another project folder

1. Choose **File > Save > Save Batch As** from the Batch View main menu in the Analysis mode.

The [Save Batch As Dialog Box](#) opens.

2. Select a storage location.

The default storage location is C:\TraceFinderData\4.0\Projects.

3. Select or create a project folder.
4. Type a name for the new batch.

If you are saving the batch to a different folder, you must give it a unique name. You cannot overwrite an existing batch in a folder.

5. Click **Save**.

When you save the batch to a different folder, the reports reflect the original project folders and the application does not save the calibration history.

❖ To move a batch to another folder

1. Choose **File > Save > Move Batch** from the Batch View main menu in the Analysis mode.

The [Save Batch As Dialog Box](#) opens.

2. Select a storage location.

The default storage location is C:\TraceFinderData\4.0\Projects.


3. Select or create a project folder.
4. Type a name for the new batch.

You must give the batch a unique name in the new subproject folder. You cannot overwrite an existing batch.

5. Click **Save**.

When you move the batch, the reports reflect the original project and subproject folders and the application does not save the calibration history.

❖ To create a new project folder

1. In the Save Batch As dialog box, select the folder for which you will create a new lower-level folder.
 - You can select the main TraceFinderData\4.0\Projects folder and create a new folder under it.
 - You can select one of the existing folders and create a lower-level folder under it.
2. Click the **Create Folder** icon, .


The application adds a new lower-level folder to the selected folder.

3. Select the new folder name and type a name for the folder.

Folder names can contain spaces and special characters, except for the following special characters: \ / : + ? " < >

Note After you add a lower-level folder, you cannot rename the parent folder.


❖ To delete project folders

1. In the Save Batch As dialog box, select the folder to delete.
2. Click the **Delete Folder** icon, .

With no confirmation prompt, the application immediately removes the selected folder.

Note This feature is not available for folders that contain lower-level project or batch folders; you must first delete the lower-level project or batch folders.

❖ To rename project folders

1. In the Save Batch As dialog box, select the folder to rename.
2. Click the **Rename Folder** icon, .

Note This feature is not available for folders that contain lower-level project or batch folders; you must first delete the lower-level project or batch folders.

3. Type a new name for the folder and press ENTER.

The application saves the new folder name.

Save Batch As Dialog Box

Use the features in the Save Batch As dialog box to save a batch to a new name or to move a batch to a different project folder.

Figure 27. Save Batch As dialog box

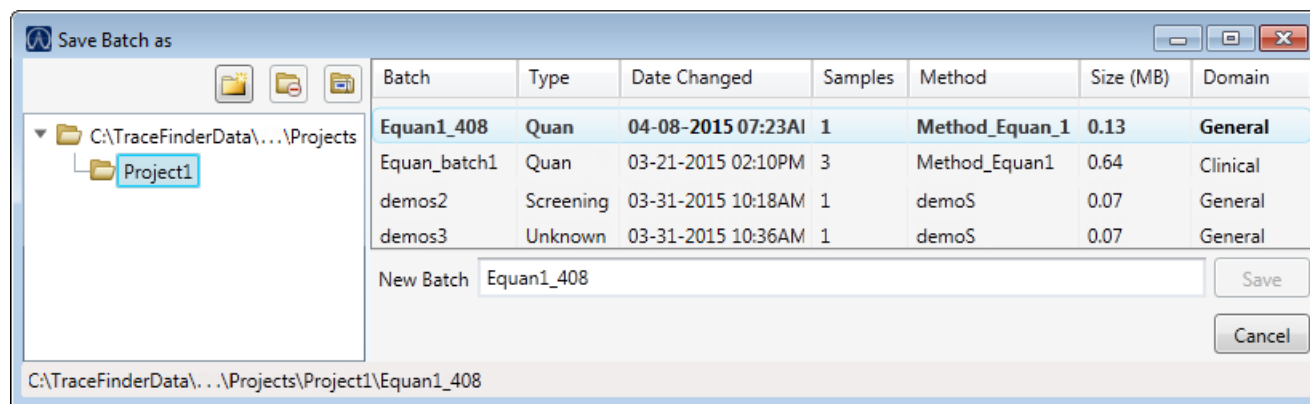


Table 25. Save Batch As dialog box parameters (Sheet 1 of 2)




Parameter	Description
 Create New Folder	Adds one of the following: <ul style="list-style-type: none"> When a drive is selected, adds a new project-level folder to the drive. When a project folder is selected, adds a subproject-level folder to the selected project. When a subproject folder is selected, adds a lower-level folder to the subproject. Or, you can right-click and choose Create Folder.
 Delete Folder	With no confirmation prompt, immediately removes the selected folder. You cannot delete a folder that contains lower-level project or batch folders; you must first delete the lower-level project or batch folders. Or, you can right-click and choose Delete.
 Rename Folder	Renames the selected folder. You cannot rename a folder that contains lower-level project or batch folders; you must first delete the lower-level project or batch folders. Or, you can right-click and choose Rename.
Batch table	
Batch	Name of batches in the selected project.
Type	Type of batch: Quan, Screening, or Unknown Only.
Date Changed	Date that the batch was last updated.
Samples	Number of samples in the batch.
Method	Name of the method used to create the batch.

Table 25. Save Batch As dialog box parameters (Sheet 2 of 2)

Parameter	Description
Size	Size of the batch in megabytes.
Domain	TraceFinder domain in which the batch was created.
New batch parameters	
New Batch	Name of the new batch to create.
	Note If the Create button is not activated, you have entered a name that is already used or you have not selected a method.
<i>Path</i>	Path to the project in the TraceFinderData\4.0\Projects folder where the batch is created.
Buttons	
Save	Saves the batch to the specified name and folder and opens the Batch View for the new batch.
Cancel	Closes the Save Batch As dialog box without saving the batch.
Shortcut menu commands	
Create Folder	<p>Adds one of the following:</p> <ul style="list-style-type: none"> • When a drive is selected, adds a new project-level folder to the drive. • When a project folder is selected, adds a subproject-level folder to the selected project. • When a subproject folder is selected, adds a lower-level folder to the subproject.
Delete Folder	<p>Immediately removes the selected folder. There is no prompt to confirm that you want to delete the selected folder.</p> <p>You cannot delete a folder that contains lower-level project or batch folders; you must first delete the lower-level project or batch folders.</p>
Rename Folder	<p>Renames the selected folder.</p> <p>You cannot rename a folder that contains lower-level project or batch folders; you must first delete the lower-level project or batch folders.</p>
Expand Child Nodes	Expands all project and subproject folders in the Project tree.
Collapse Child Nodes	Collapses all project and subproject folders in the Project tree.

Auto Samples Page for Quantitation Batches

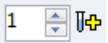
The Auto Samples page identifies the Solvent or Negative samples to use for any Auto Sample or Auto Sample and Reinject failure actions as specified on the Intelligent Sequencing page of the method. Refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

Each sample type that you specify for a failure action on the Intelligent Sequencing page must be defined in the samples list on the Auto Samples page.

❖ To add an auto sample type

1. Click **Auto Samples** in the Batch View navigation pane.

The [Auto Samples page](#) opens.

2. Right-click and choose **Add Auto Sample**, or in the toolbar click the **Add New Auto Sample** icon, .

The application adds a Solvent sample to the sample list.

You can add, insert, or remove samples from this list as you would any sample list. See [Samples Page for Quantitation Batches](#).

3. To change the sample type to a Negative, click the Sample Type column and select Negative from the list.
4. In the Injection Volume column for the sample, type a volume.

The minimum injection volume value allowed is 0.1 µL; the maximum injection volume value allowed is 5000 µL.

5. In the Number of Injections column, type the number of injections available in the designated Solvent or Negative vial.

After auto sample injections have occurred, you can return to this page to view the number of Injections Used in each vial.

6. In the Vial Position column, type the vial position for the Solvent or Negative sample.

Auto Samples Page

Use the Auto Samples page to specify the sample types to use with Intelligent Sequencing.

Figure 28. Auto Samples page

	Sample Type	Injection Volume	Injections Used	Number of Injections	Vial Position
	Solvent	1.0	0	1	10
	Matrix Blank	1.0	0	10	11
	Matrix Blank	1.0	0	10	12

Table 26. Auto Samples page parameters

Column	Description
Sample Type	The sample type for the auto sample injection as specified on the Intelligent Sequencing page—either Solvent or Negative. Default: Solvent
Injection Volume	The injection volume used for the sample acquisition as specified on the Samples page. Valid range: 0.1 through 5000 µL
Injections Used	The number of times a vial has been used. The count is cumulative across all batches.
Number of Injections	The number of injections available in the designated Solvent or Negative vial.
Vial Position	Vial position for this sample type as specified on the Samples page.

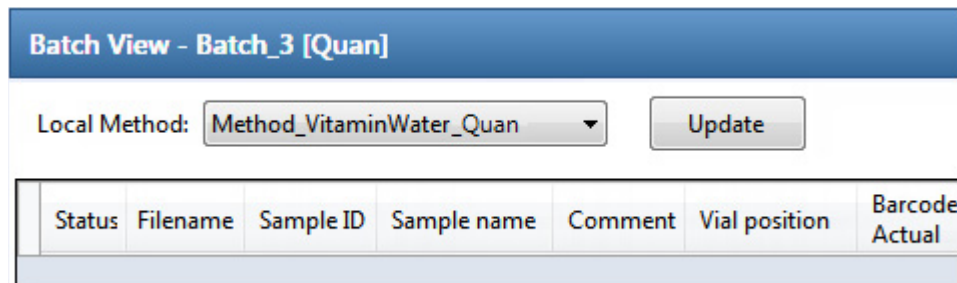
Reference Sample Page for Quantitation Batches

The Reference Sample page displays the reference samples that you selected for this batch.

❖ To specify a chromatogram reference sample

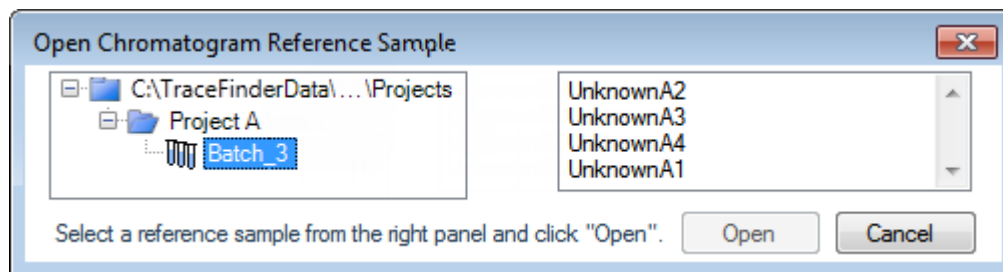
1. In the Batch View, click **Reference Sample**.

An empty reference sample table opens.



2. Click the **Add Reference Sample** icon, , in the toolbar, or right-click and choose **Add Reference Sample**.

The Open Chromatogram Reference Sample dialog box opens.



Note If you are using a new method, no reference samples appear here. You must first process a batch using the current method to see the reference samples in this list.

3. Select a project from the list of projects.
4. Select a subproject from the list of subprojects.
5. Select a batch from the list of batches.

The application displays only batches that were created using the current master method.

6. Select a sample from the list of processed samples.

The application displays all the processed samples in the selected batch. Before using a sample as a reference sample, you must have processed the sample with the current master method.

7. Click **Open**.

Threshold Samples Page for Quantitation Batches

For each group in a batch, you can specify a sample in the group as the threshold sample to use in the Comparative View.

❖ To specify a threshold sample

1. In the Batch View, click **Threshold Samples**.
2. Click the Sample list for each group and select a sample in the group to be the threshold sample.

	Group	Sample
▶	groupb	Benzo26473 ▼
		Benzo26473
		Benzo25557
		Benzo26154

The Comparative View uses the threshold method and amount you specified in the method, the group you created on the Samples page, and the threshold sample you selected on this page to define the threshold guide that it displays in the sample peak plots.

To specify the method for creating a threshold guide, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

For information about creating groups, see [Groups](#).

For information about using the threshold guide in the Comparative View, see [Comparative View for Quantitation Batches](#).

Working in Data Review for Quantitation Batches

In the Data Review view, you can view the data generated by the master method. Use Data Review to verify the data for a compound before you generate reports.

Follow these procedures:

- [To open the Data Review view](#)
- [To move, dock, or float Data Review panes](#)
- [To restore the default layout](#)

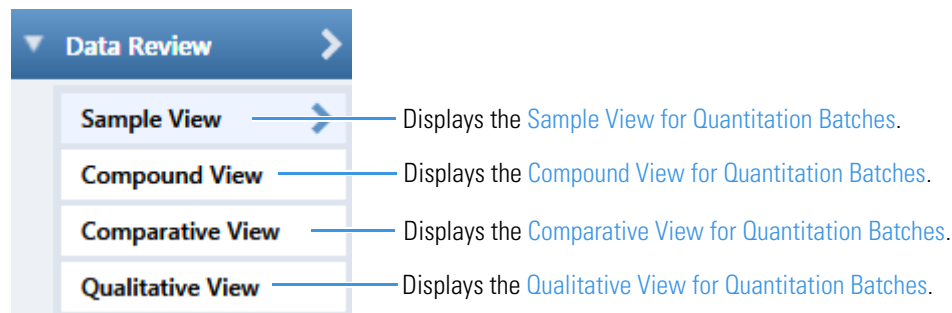
❖ To open the Data Review view

1. Click **Analysis** in the navigation pane.

The Analysis navigation pane opens.

2. Click **Data Review**.

The Data Review navigation pane opens.



Choose from a Sample View, Compound View, Comparative View, or Qualitative View to analyze the data generated by the master method.

In addition to these views, see the following topics:

- [Compound Details for Quantitation Batches](#)
- [Column Parameters for Compound Results and Sample Results](#)

❖ To move, dock, or float Data Review panes

Follow the instructions in [Appendix C, “Moving Data Review Panes.”](#)

❖ To restore the default layout

Choose **View > Restore Default Layout**.

The Sample View, Compound View, Comparative View, and Qualitative View have their own defaults for which panes are displayed and in which location.

Sample View for Quantitation Batches

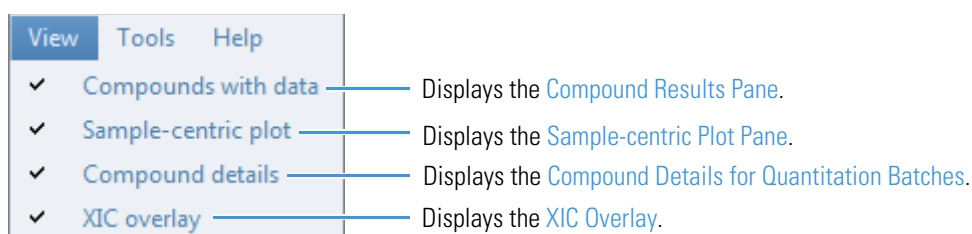
The Sample View displays a list of all samples in the current batch and the compound results for all compounds in the method.

❖ To display or hide a pane on the Sample View page

From the View menu, choose to display or hide any of the following panes.

Note The Samples pane is required for the Sample View display. You cannot hide the Samples pane.

A check mark indicates a displayed pane.



The Sample View display includes the following features:

- [Samples Pane](#)
- [Compound Results Pane](#)
- [Sample-centric Plot Pane](#)
- [XIC Overlay](#)

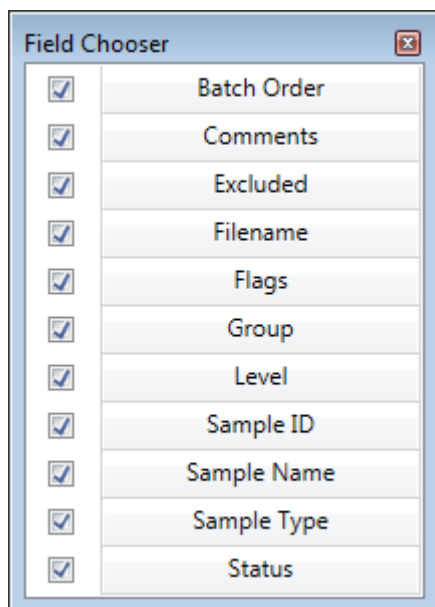
Samples Pane

Use the Samples pane in the Sample View to select a specific sample. The associated [Compound Results Pane](#) displays all compounds in the method and flags any compound with errors in the selected sample.

❖ To hide or display columns in the Samples pane

1. Click the **Field Chooser** icon, , in the upper left corner of the pane.

The Field Chooser displays all available columns of data for the Samples pane.



Note The Field Chooser also lists any custom columns that you defined in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.

2. Select the check box for each column that you want to display, or clear the check box for each column that you want to hide.

The application immediately displays or hides the column in the Samples pane.


3. When you are finished modifying the column display, click  to close the Field Chooser.

Figure 29. Samples pane

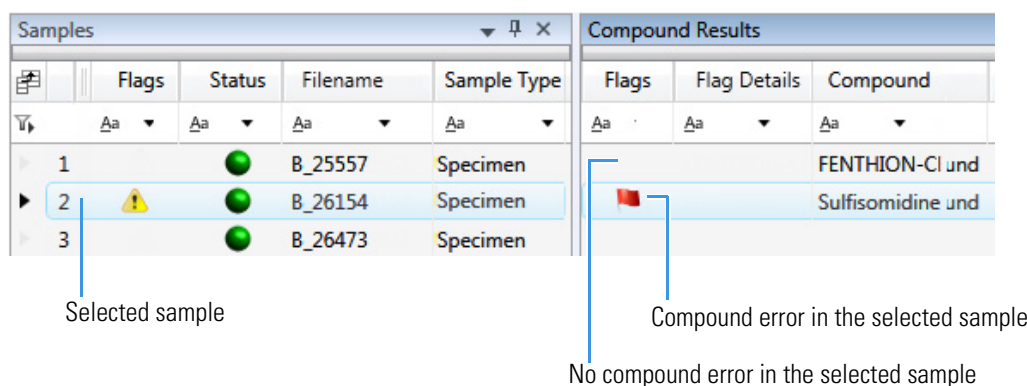




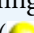





Table 27. Samples pane columns

Column	Description
Batch Order	Sequentially numbers the samples.
Comments	User-defined comments for the sample.
Excluded	Turns a compound on or off in the calibration curve in the Compound Details pane.
Filename	A user-defined name that identifies a sample.
Flags	Caution flag displayed when a compound in the sample has an error. See Caution Flags .
Group	Threshold group to which a sample belongs. Samples can be viewed by group in the Comparative View of Data Review.
Level	The level defined for a calibration sample or quality control sample.
Sample ID	A user-defined, alphanumeric string that identifies a sample.
Sample Name	A user-defined name that identifies a sample.
Sample Type	Defines how the application processes the sample data. Each sample is classified as one of the following sample types: Specimen, QC, Solvent, Calibrator, Hydrolysis, Unextracted, or Negative.
Status	 Sample is not acquired.  Sample is acquired but not processed.  Sample is acquired and processed.  Sample is currently acquiring.
Note When you include unknown screening features in the quantitation method and you choose to process with only the quantitation criteria, the Sample View shows the Status for the samples as acquired and processed () , and the Unknown Screening View shows the Status for the samples as acquired but not processed ().	
Note When you include unknown screening features in the quantitation method and you choose to process with only the unknown screening criteria, both the Sample View and the Unknown Screening View show the Status for the samples as acquired and processed ().	

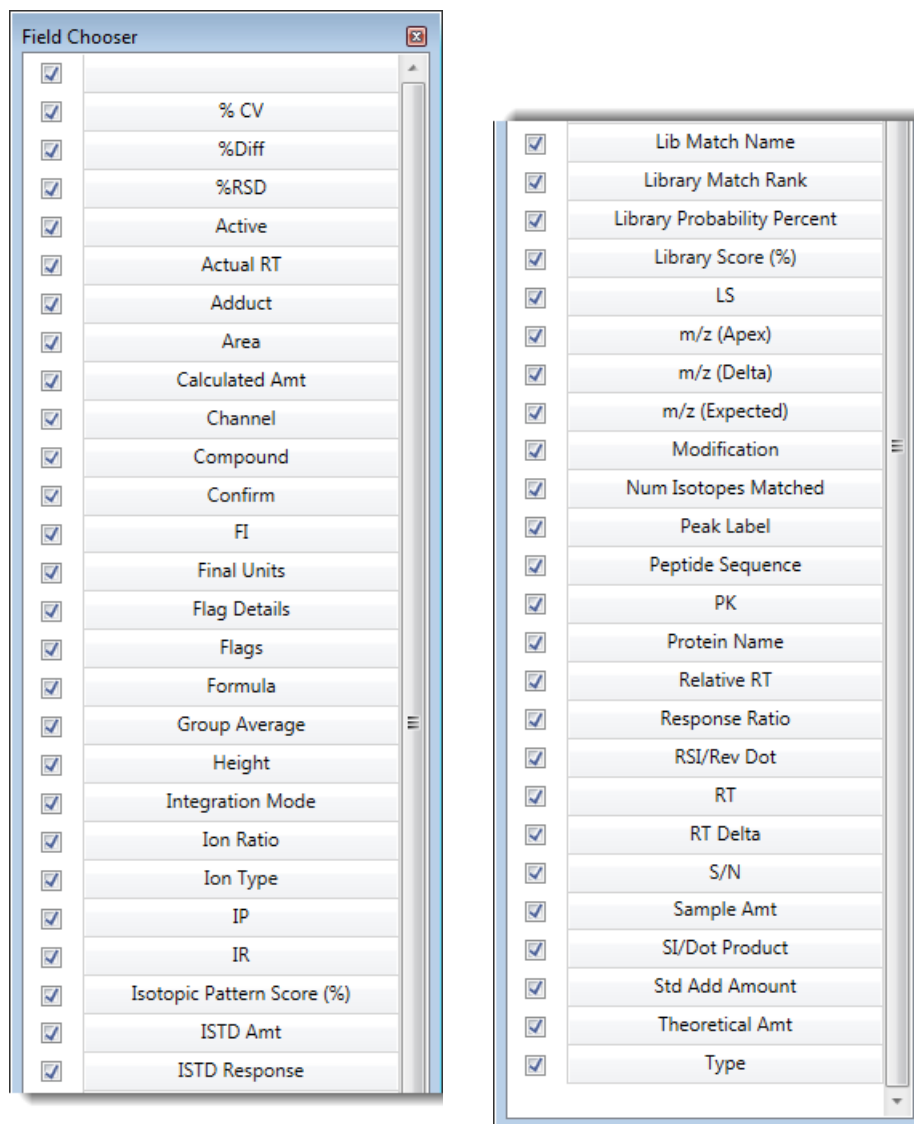
Compound Results Pane

Use the Compound Results pane in the Sample View to select a specific compound in the selected sample. The associated [Sample-centric Plot Pane](#) highlights the selected compound.

❖ To hide or display columns in the Compound Results pane


1. Click the **Field Chooser** icon, , in the upper left corner of the pane.

The Field Chooser displays all available columns of data for the Compound Results pane.



2. Select the check box for each column that you want to display, or clear the check box for each column that you want to hide.

The application displays or hides the column in the Compound Results pane.

3. When you are finished modifying the column display, click  to close the Field Chooser.

4 Using the Analysis Mode for Quantitation Batches

Working in Data Review for Quantitation Batches

Figure 30. Compound Results pane

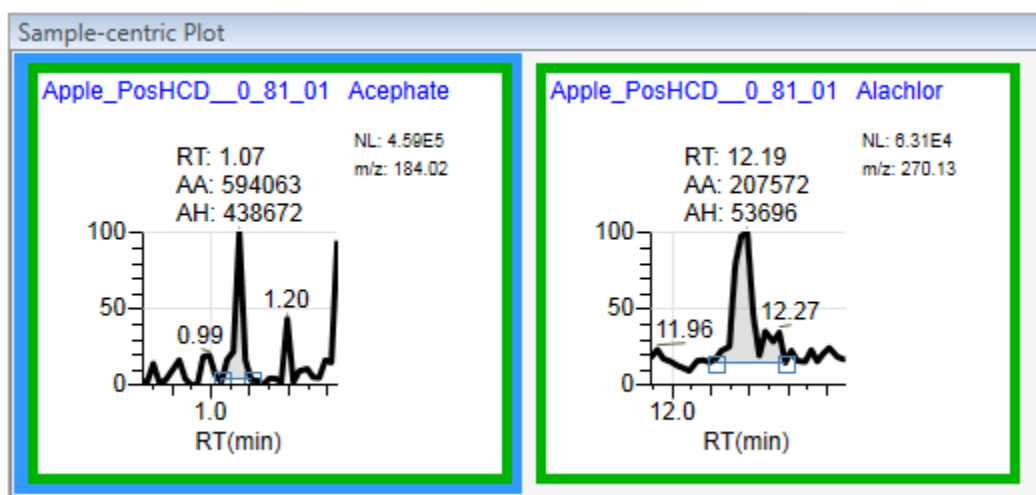
Compound Results							
	Flags	Flag Details	Peak Label	Compound	Type	Height	Area
	Aa	Aa	Aa	Aa	Aa	Aa	Aa
1	<input checked="" type="checkbox"/>			Acephate	Target Compound	2537075	1398474
1	<input checked="" type="checkbox"/>		T1	Acephate	Target Peak	396175	376065
2	<input checked="" type="checkbox"/>		T1F1	Acephate	Fragment	N/F	N/A
3	<input checked="" type="checkbox"/>		T1F2	Acephate	Fragment	N/F	N/A
4	<input checked="" type="checkbox"/>		T1F3	Acephate	Fragment	N/F	N/A
5	<input checked="" type="checkbox"/>		T2	Acephate	Target Peak	2140900	102240
6	<input checked="" type="checkbox"/>		T2F1	Acephate	Fragment	N/F	N/A
7	<input checked="" type="checkbox"/>		T2F2	Acephate	Fragment	N/F	N/A
8	<input checked="" type="checkbox"/>		T2F3	Acephate	Fragment	N/F	N/A

For descriptions of the column parameters in the Compound Results pane, see [Common Column Parameters](#).

Sample-centric Plot Pane

The Sample-centric Plot pane in the Sample View displays the chromatogram, retention time, area, height, and signal-to-noise ratio for each compound in the Compound Results pane. The application highlights the chromatogram for the compound that is currently selected in the Compound Results pane.

You can use the Sample-centric Plot pane to view an overlay of the quantitation and confirming peaks for each sample in the batch.



Follow these procedures:

- To select a sample to view
- To display details for a compound
- To change the number of rows and columns in the display

- To change the target and confirming peaks that are displayed
- To zoom in on a peak
- To copy a plot to another format

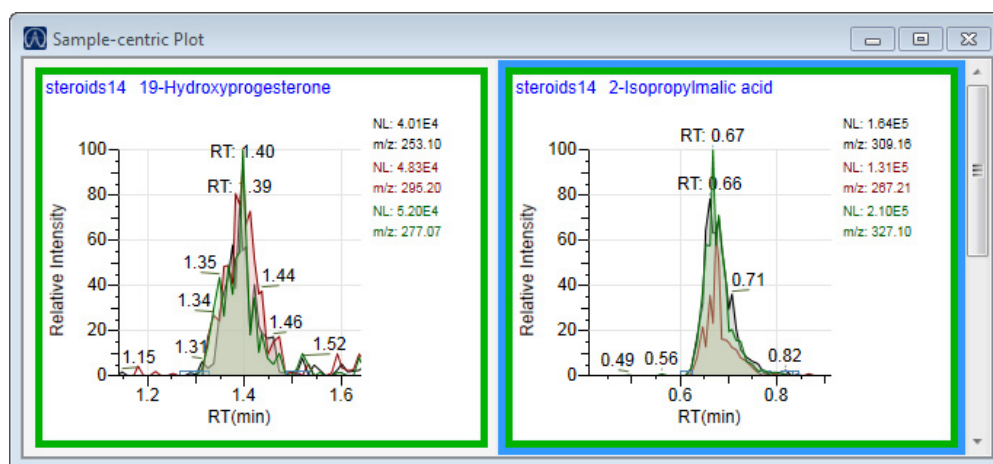
❖ **To select a sample to view**

1. In the Samples pane, select the sample that you want to view.

Samples				Compound Results	
	Bat	Status	Filename	Flags	Compound
1	1	●	steroids02	1	19-Hydroxyprogesterone
2	2	●	steroids04	2	2-Isopropylmalic acid
3	3	●	steroids14	3	Glycitein
4	4	●	steroids15	4	Progesterone

The Sample-centric Plot pane displays the compound chromatograms for all compounds in the selected sample. A blue border indicates the compound in the sample that is currently selected in the Compound Results pane.

Figure 31. Sample-centric Plot pane



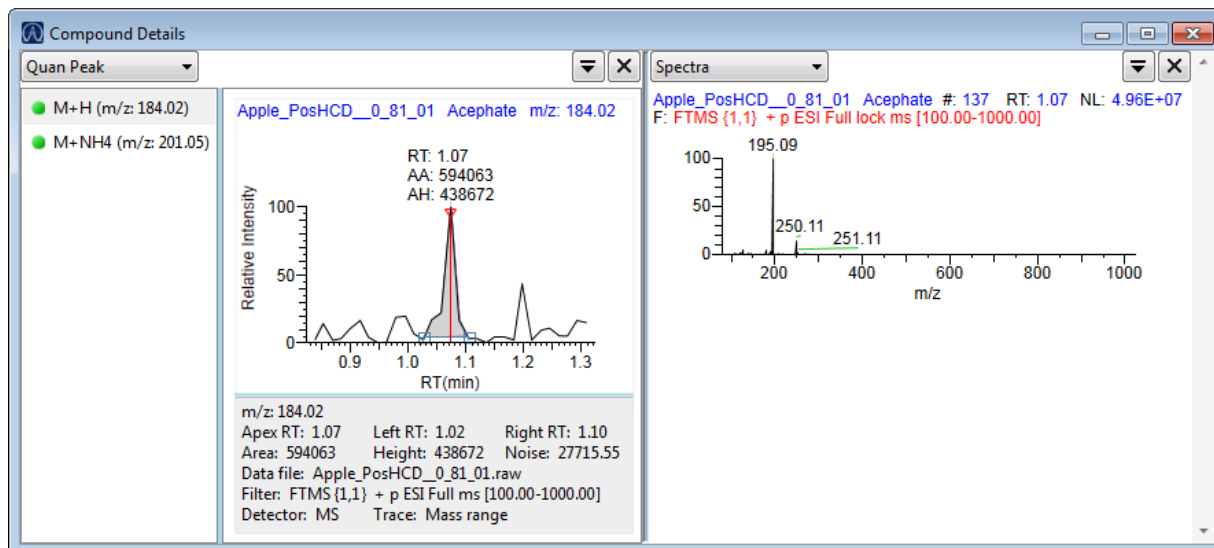
4 Using the Analysis Mode for Quantitation Batches

Working in Data Review for Quantitation Batches

❖ To display details for a compound

Double-click the chromatogram in the Sample-centric Plot pane.

The Compound Details pane opens.



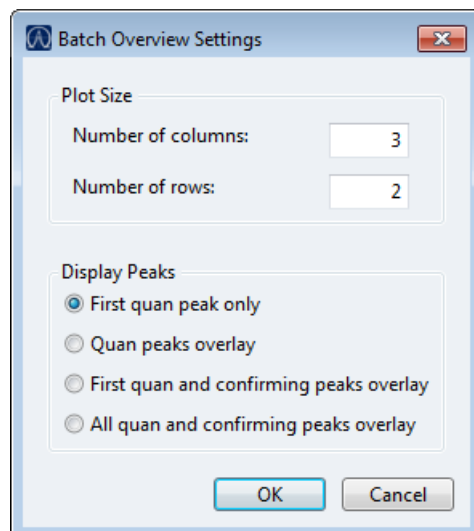
The Compound Details pane displays information about the quantitative peak, calibration curve, confirming ion, internal standard, reference peak, ion overlay, and spectra for the compound.

For a detailed description of the available information in the Compound Details pane, see [Compound Details for Quantitation Batches](#).

❖ To change the number of rows and columns in the display

1. Right-click and choose **Display Settings**.

The Batch Overview Settings dialog box opens.



- To change the number of rows or columns to fit in the Sample-centric Plot pane, type new values in the Number of Columns or Number of Rows boxes.

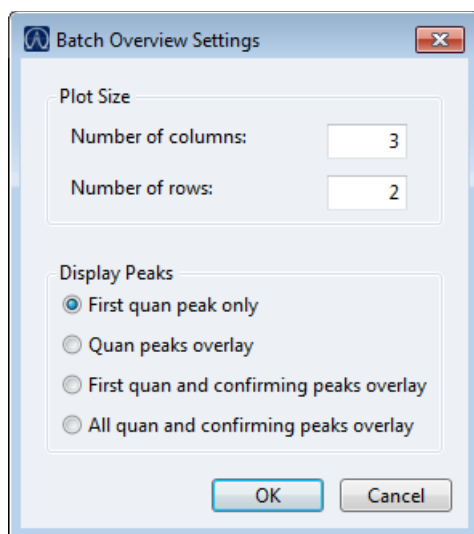
These values do not change the number of rows (compounds) and columns (samples) that are available in the Sample-centric Plot pane. They determine how many rows and columns you want to view at one time in the display. The default is two rows and three columns. Use the scroll bar to view all compound chromatograms.

- Click **OK**.

❖ **To change the target and confirming peaks that are displayed**

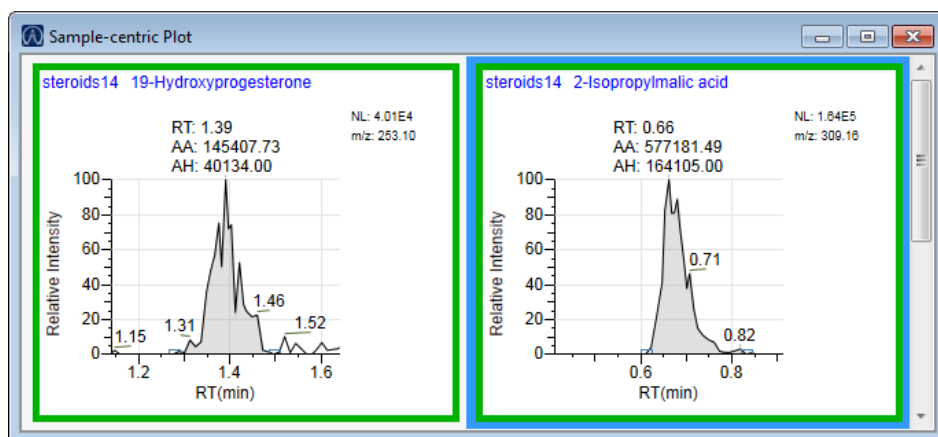
- Right-click and choose **Display Settings**.

The Batch Overview Settings dialog box opens.



- Select any of the following display options:

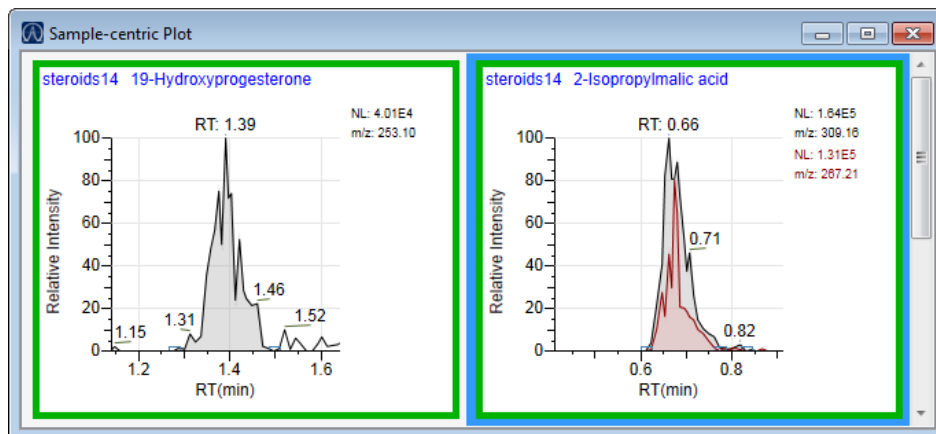
- First Quan Peak Only



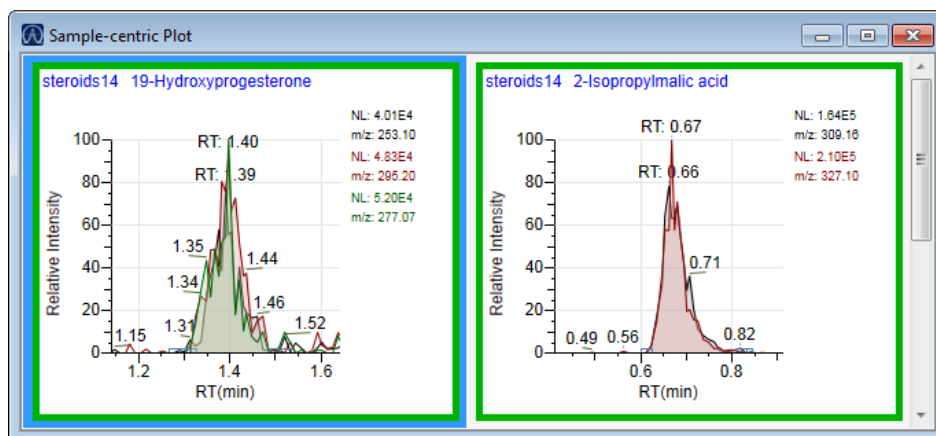
4 Using the Analysis Mode for Quantitation Batches

Working in Data Review for Quantitation Batches

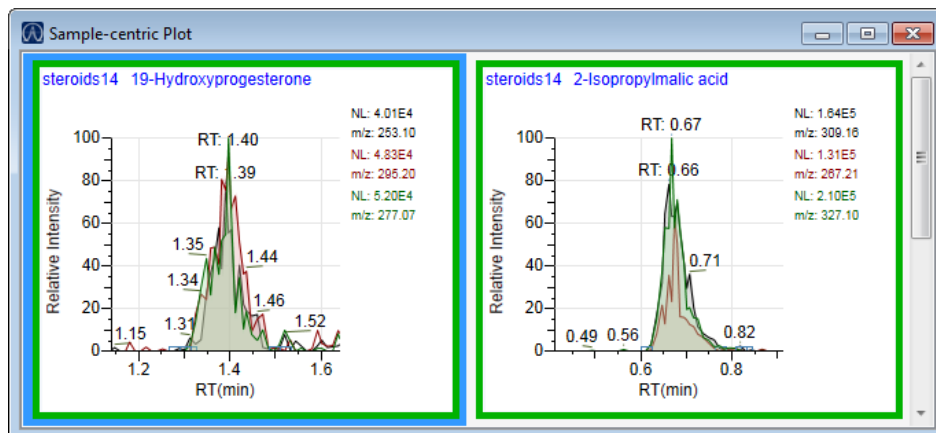
- Quan Peaks Overlay



- First Quan and Confirming Peaks Overlay



- All Quan and Confirming Peaks Overlay



3. Click **OK**.

❖ **To zoom in on a peak**

1. In any of the chromatogram cells, drag your cursor to delineate a rectangle around the peak.

The delineated area expands to fill the chromatogram area.

2. To restore the default view, right-click the chromatogram plot and choose **Reset Scaling**.

❖ **To copy a plot to another format**

1. Right-click the plot that you want to copy and choose **Copy to Clipboard**.

The application copies the plot graphic to the Clipboard.

2. In the source format (for example, an Excel spreadsheet or email window), right-click and choose **Paste**.

The application pastes the graphic into the new format.

XIC Overlay

Use the XIC Overlay pane to view specific groups of peaks. You can choose to view all peaks, selected peaks only, or the top 20 most intense peaks (by area). The XIC Overlay plot is a collection of overlaid, extracted m/z ion plots that use a different color for each peak.

Follow these procedures:

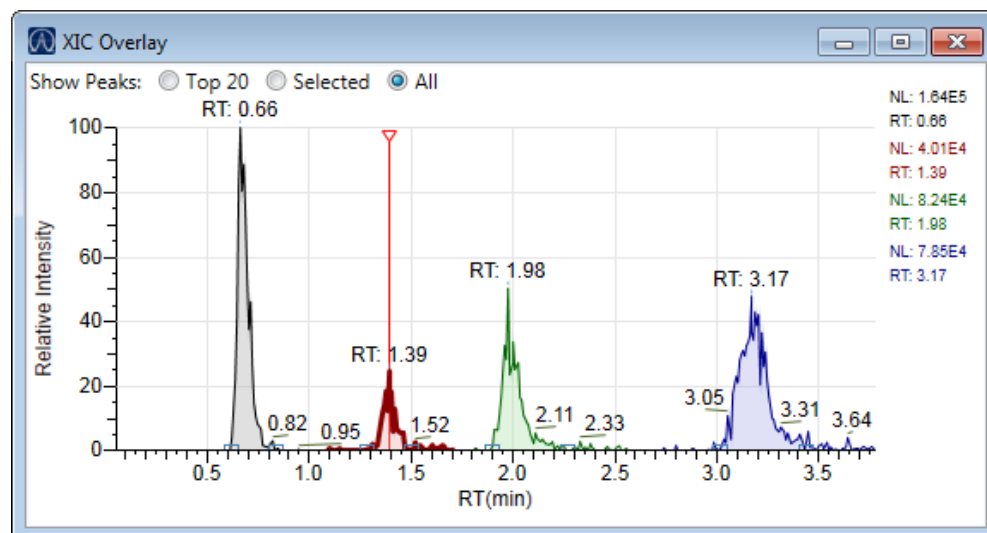
- To display all peaks in a sample
- To display only selected peaks in a sample
- To display the top 20 peaks in a sample
- To focus on a specific peak

❖ **To display all peaks in a sample**

1. In the Samples pane, select the sample whose peaks you want to view.
2. In the XIC Overlay pane, select the **All** check box.

The application displays all the peaks in the selected sample.

Figure 32. XIC Overlay pane showing all peaks



Note If there are more than 1000 peaks in the sample, performance might be slow.

❖ To display only selected peaks in a sample

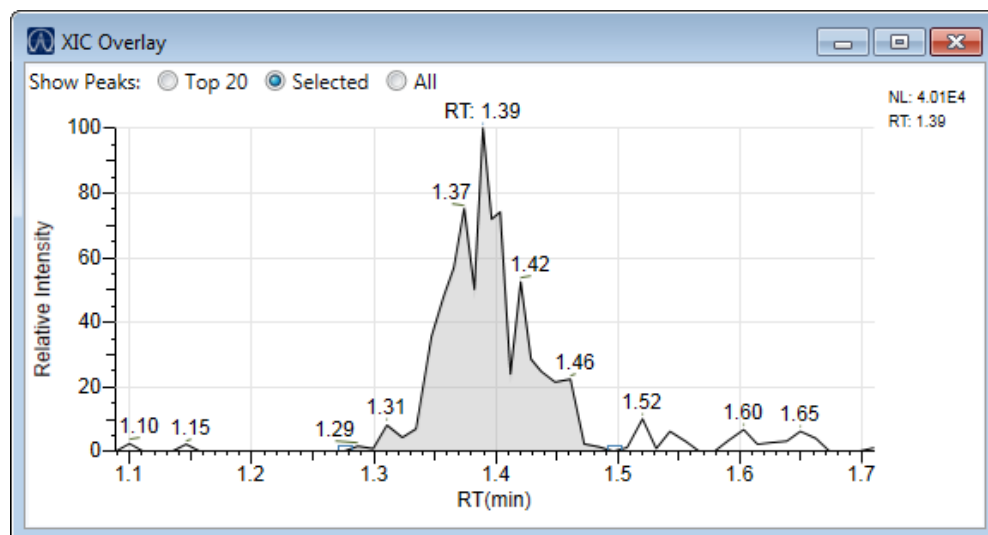
1. In the Samples pane, select the sample whose peaks you want to view.
2. In the Compound Results pane, select the check box for each peak that you want to view.

Compound Results				
			Compound	Type
1	<input checked="" type="checkbox"/>		19-Hydroxyprogesterone	Target Compound
1	<input checked="" type="checkbox"/>		19-Hydroxyprogesterone	Target Peak
2	<input checked="" type="checkbox"/>		19-Hydroxyprogesterone	Confirming Peak
3	<input checked="" type="checkbox"/>		19-Hydroxyprogesterone	Confirming Peak

3. In the XIC Overlay pane, choose the **Selected** option.

The application displays only the selected peaks.

Figure 33. XIC Overlay pane showing selected peaks

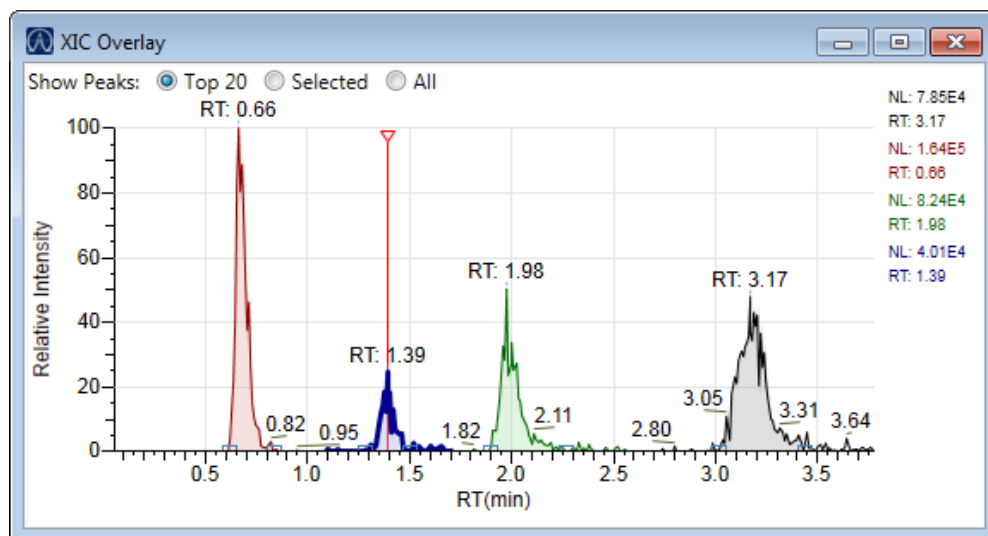


❖ **To display the top 20 peaks in a sample**

1. In the Samples pane, select the sample whose peaks you want to view.
2. In the XIC Overlay pane, select the **Top 20** option.

The application displays the top 20 peaks in the selected sample.

Figure 34. XIC Overlay pane showing the top 20 peaks



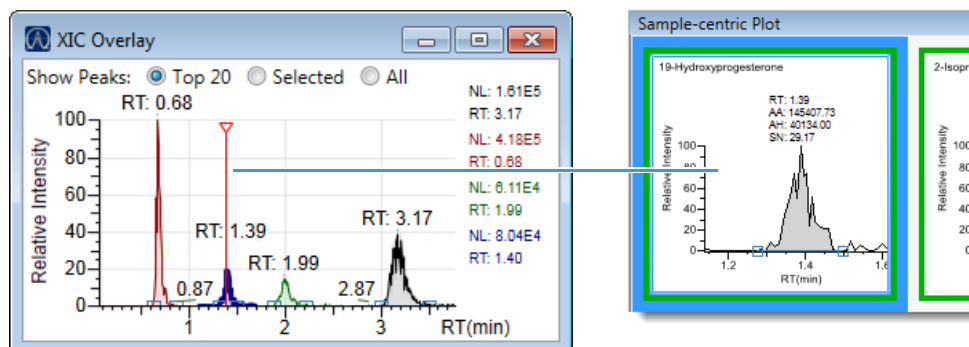
4 Using the Analysis Mode for Quantitation Batches

Working in Data Review for Quantitation Batches

❖ To focus on a specific peak

In the XIC Overlay pane, click any displayed peak.

The application indicates the selected peak with a red, vertical marker and updates the focus in the other Data Review panes with data for that peak.



Caution Flags


In the Sample View, the application displays caution flags in both the Samples pane and the Compound Results pane.

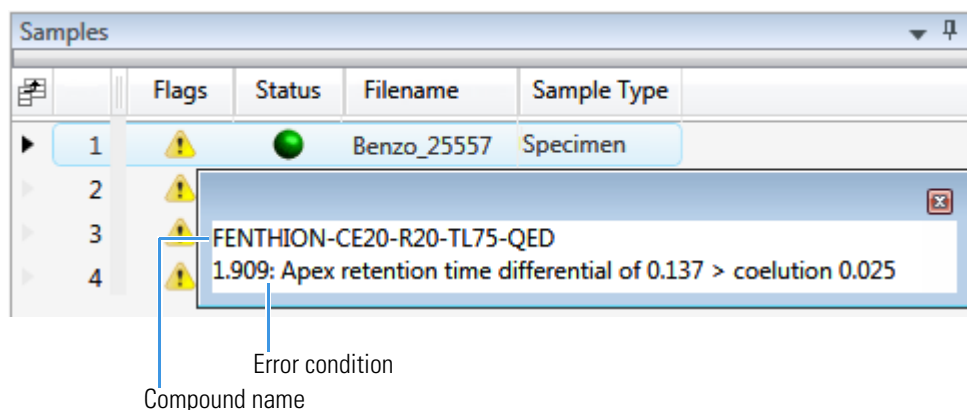
See the following topics:

- [Flags in the Samples Pane](#)
- [Flags in the Compound Results Pane](#)
- [Error Indicators in the Sample-centric Plot Pane](#)

Flags in the Samples Pane

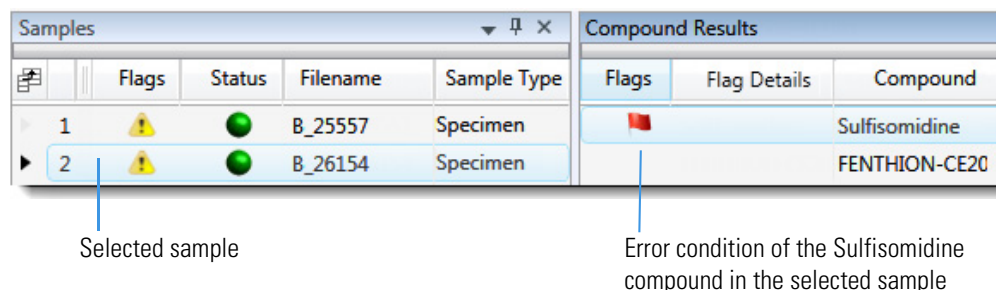
The Flags column in the Samples pane displays a caution flag if any compound in the sample is not in compliance with the method criteria.


Click the caution flag icon, , to display the details. Information in the pop-up box shows the compound that is in error and describes the exact error condition.

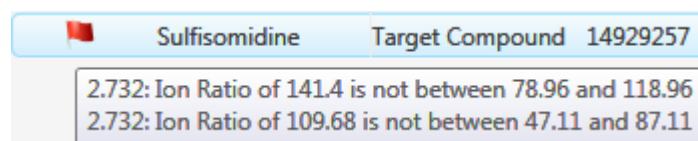


Flags in the Compound Results Pane







The Flags column in the Compound Results pane displays a flag if the compound in the selected sample is not in compliance with the method criteria.



Hold your cursor over the flag icon, , to display details for the compound in the selected sample.



Colored flags in the Compound Results pane have the following meaning:

-  Red for compounds that have violated (or are activated by) any of the values set in the method. Refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.
-  Red for compounds that are outside the specified ion ratio range. See [Ion ratio failure flag](#).
-  Orange for compounds that are below the limit of quantitation (LOQ), below the limit of detection (LOD), or between the LOD and LOQ values specified in the method. For descriptions of these limits, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.
-  Green for “found” compounds that are over the limit of reporting (LOR) amount specified in the method. For a description of the LOR limit, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.
-  Yellow for compounds that are equal to or below the LOR amount specified in the method.
-  Yellow for compounds that are not found in Calibrator or QC sample types. The Compound Results pane does not flag compounds that are not found in Specimen sample types.

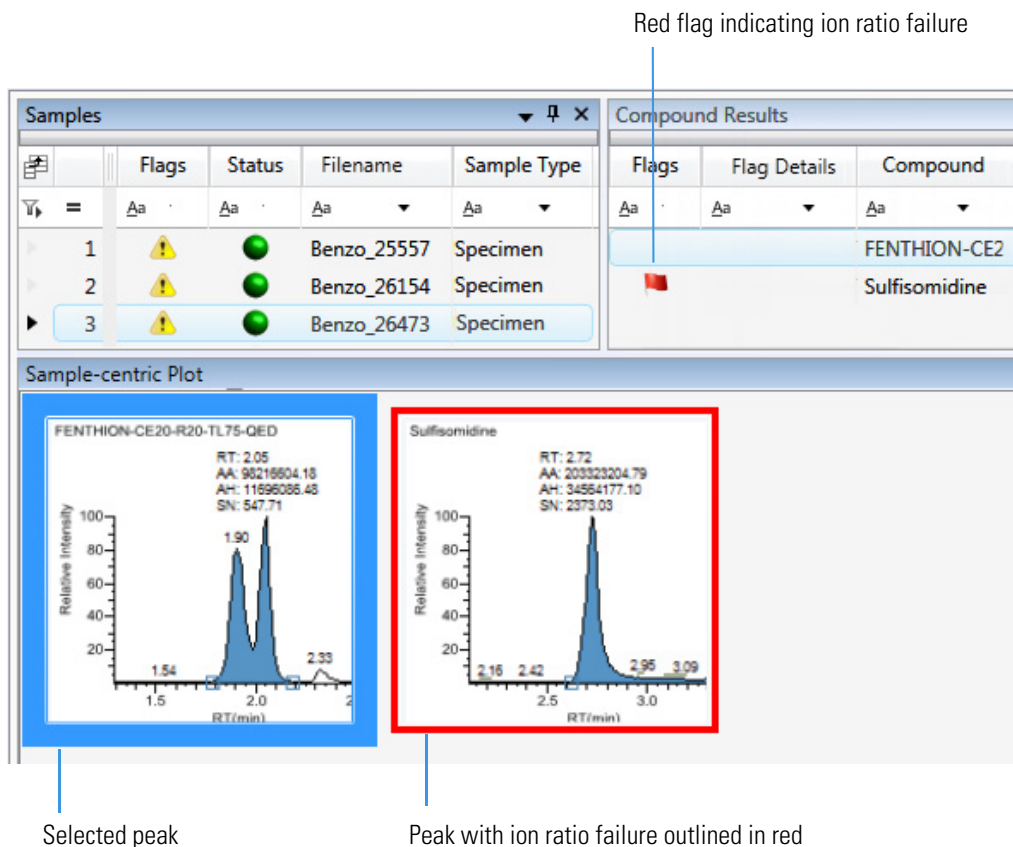
No flag for compounds that have no errors or where no report options are selected.

Note These criteria for flag states do not apply to Negative sample types when the compound is an internal standard.

Error Indicators in the Sample-centric Plot Pane

In the Sample-centric Plot pane, peak plots are outlined with the color of their associated error flag. In the following example, the FENTHION peak plot is highlighted in blue to indicate that FENTHION is the selected compound, and the Sulfisomidine peak plot is outlined in red to indicate that the Sulfisomidine compound in the selected sample is outside the specified ion ratio range.

Figure 35. Ion ratio failure flag



Compound View for Quantitation Batches

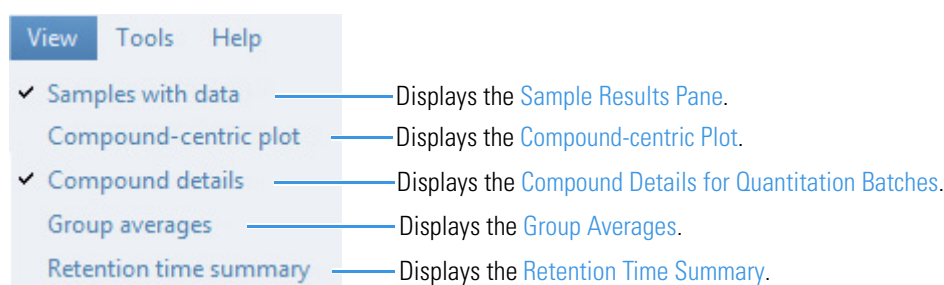
The Compound View displays a list of all compounds available in the method, all samples in the current batch, and the peak plots for all found compounds.

❖ To display or hide a pane on the Compound View page

From the View menu, choose to display or hide any of the following panes.

Note The Compounds pane is required for the Compound View display. You cannot hide the Compounds pane.

A check mark indicates a displayed pane.




The Compound View includes the following features:

- [Compounds Pane](#)
- [Sample Results Pane](#)
- [Compound-centric Plot](#)
- [Group Averages](#)
- [Retention Time Summary](#)
- [Caution Flags](#)

Compounds Pane

Use the Compounds pane in the Compound View to select a specific compound. The [Sample Results Pane](#) displays all samples in the batch and flags any sample that contains errors associated with the selected compound.

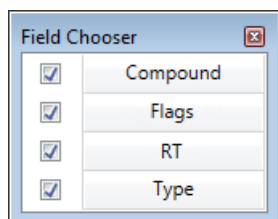
❖ To hide or display columns in the Compounds pane

1. Click the **Field Chooser** icon, , in the upper left corner of the pane.

4 Using the Analysis Mode for Quantitation Batches

Working in Data Review for Quantitation Batches

The Field Chooser displays all available columns of data for the Compounds pane.



2. Select the check box for each column that you want to display, or clear the check box for each column that you want to hide.

The application immediately displays or hides the column in the Compounds pane.


3. When you are finished modifying the column display, click  to close the Field Chooser.

Figure 36. Compounds pane

Compounds				Sample Results					
Flags	Compound	RT	Type		Batch	Flags	Flag Details	Status	Filename
1	19-Hydroxyprogesterone	1.39	Target Compound	+	1	1	ICPF	●	steroids02
2	2-Isopropylmalic acid	0.68,0.68	Target Compound	+	2	2		●	steroids04
3	Glycitein	1.98	Target Compound	+	3	3	ICPF	●	steroids14
4	Progesterone	3.15	Target Compound	+	4	4		●	steroids15

Selected compound

Error associated with the selected compound in this sample

No error associated with the selected compound in this sample


Table 28. Compounds pane columns

Column	Description
Flags	Caution flag displayed when a compound has an error in any of the samples.
Compound	Compound names as identified in the library. If there is no library selected in the method template, the compound name is identified as <i>peak@RT</i> .
RT	Expected retention time for the compound.
Type	Specified compound type: Target Compound or Internal Standard.

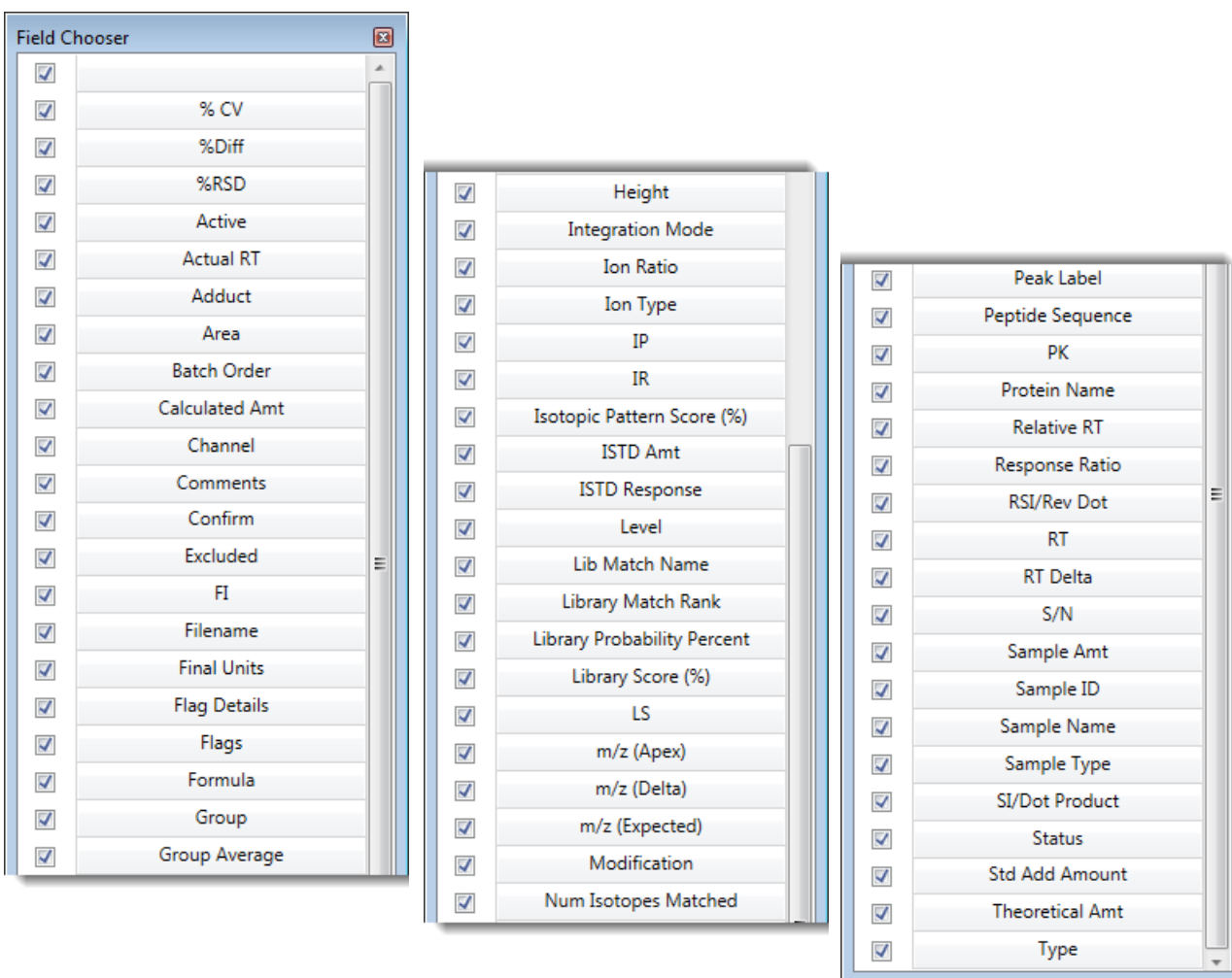
Sample Results Pane

Use the [Sample Results pane](#) in the Compound View to select a specific compound in a specific sample. The Compound-centric Plot pane highlights the selected sample/compound and displays the following information about the compound: chromatogram, retention time, area, height, and signal-to-noise ratio.

❖ To hide or display columns in the Sample Results pane


1. Click the **Field Chooser** icon, , in the upper left corner of the pane.

The Field Chooser displays all default columns of data for the Sample Results pane.



2. Select the check box for each column that you want to display, or clear the check box for each column that you want to hide.

The application immediately displays or hides the column in the Sample Results pane.

3. When you are finished modifying the column display, click  to close the Field Chooser.

4 Using the Analysis Mode for Quantitation Batches

Working in Data Review for Quantitation Batches

Figure 37. Sample Results pane

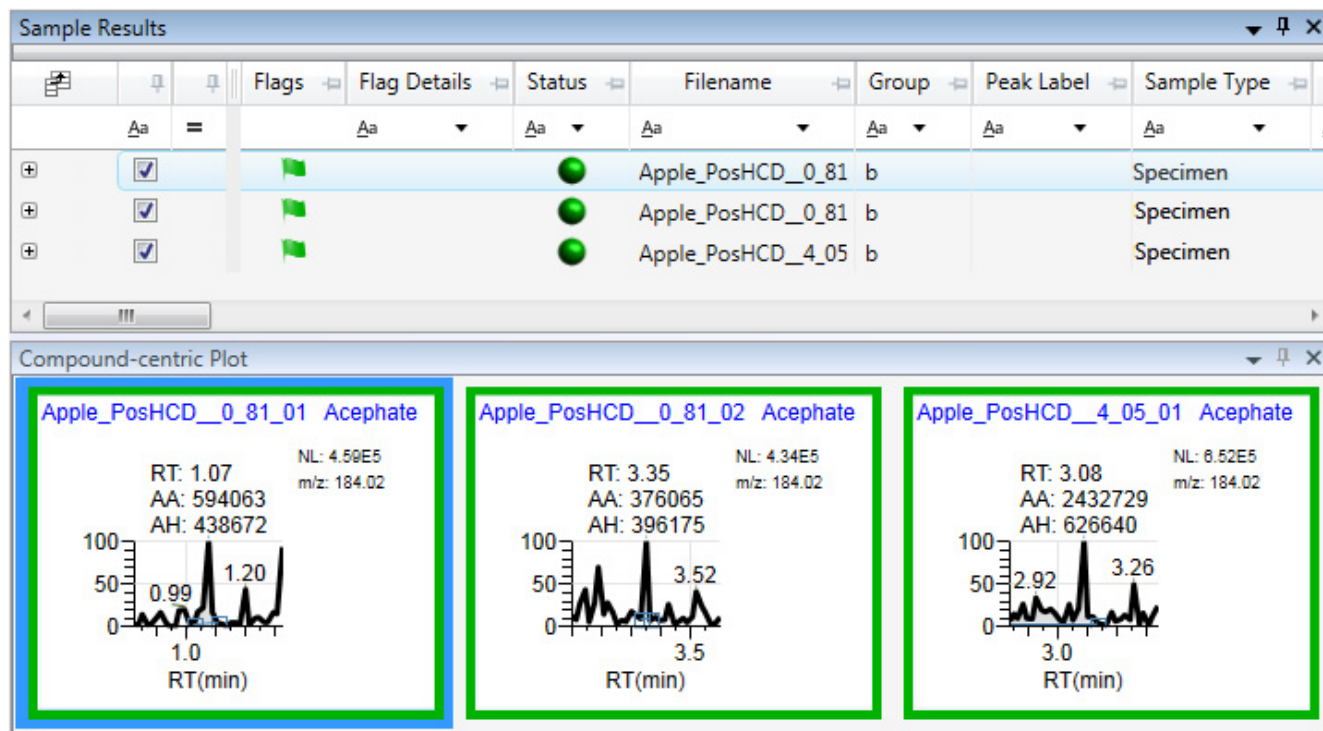


Table 29. Sample Results pane columns (Sheet 1 of 2)

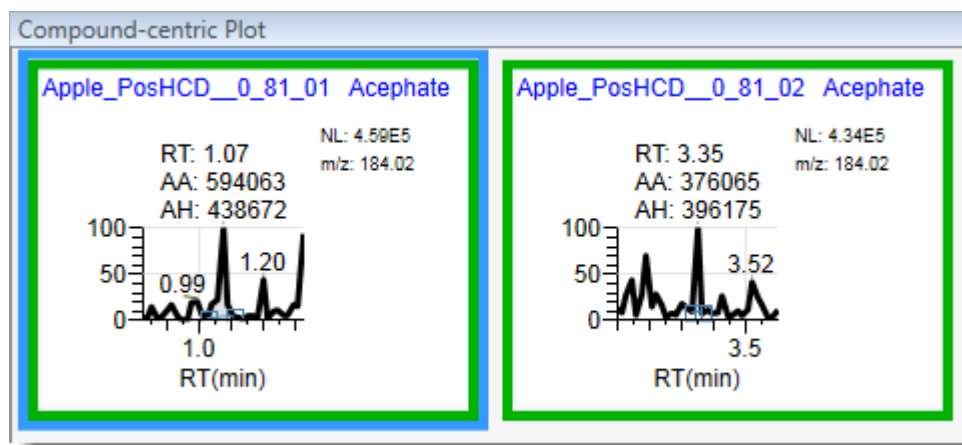
Column	Description
<input checked="" type="checkbox"/>	Displays the selected samples in the other Data Review panes.
Flags	Displays a caution flag when a compound within the sample has an error.
Flag Details	Indicates the type of error: <ul style="list-style-type: none"> • I: Confirming ion coelution failure or Ion ratio failure • A: Amount error • B: Matrix blank error • H: Peak not found

Table 29. Sample Results pane columns (Sheet 2 of 2)

Column	Description
Status	<p>● Sample is not acquired.</p> <p>● Sample is acquired but not processed.</p> <p>● Sample is acquired and processed.</p> <p>● Sample is currently acquiring.</p> <p>Note When you include unknown screening features in the quantitation method and you choose to process with only the quantitation criteria, the Sample View shows the Status for the samples as acquired and processed (●), and the Unknown Screening View shows the Status for the samples as acquired but not processed (●).</p> <p>Note When you include unknown screening features in the quantitation method and you choose to process with only the unknown screening criteria, both the Sample View and the Unknown Screening View show the Status for the samples as acquired and processed (●).</p>
Filename	Specifies a user-defined name that identifies a sample.
Group	Specifies the threshold group to which a sample belongs. Samples can be viewed by group in the Comparative View of Data Review.
Sample Type	Defines how the application processes the sample data. Each sample is classified as one of the following sample types: Specimen, QC, Solvent, Calibrator, Hydrolysis, Unextracted, or Negative.
The remainder of the columns in the Sample Results pane are common to both the Sample View and the Compound View displays. See Common Column Parameters .	

Compound-centric Plot Pane

The Compound-centric Plot pane in the Compound View displays the compound chromatogram, retention time, area, height, and signal-to-noise ratio associated with the selected compound in each of the samples in the batch. The application highlights the compound chromatogram for the sample that is currently selected in the Sample Results pane.



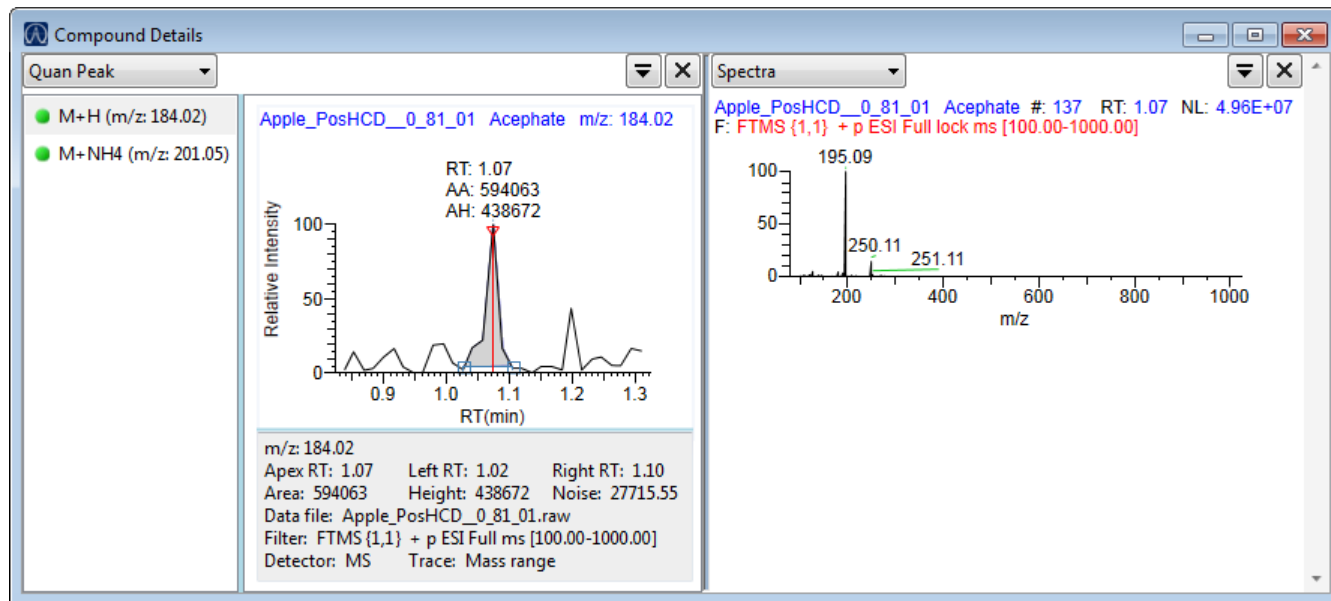
4 Using the Analysis Mode for Quantitation Batches

Working in Data Review for Quantitation Batches

❖ To display details for a compound

Double-click the chromatogram in the Compound-centric Plot pane.

The application adds the Compound Details pane to the window.



The Compound Details pane displays information about the quantitative peak, calibration curve, confirming ion, internal standard, reference peak, ion overlay, and spectra for the compound.

For details about the available information in the Compound Details pane, see [Compound Details for Quantitation Batches](#).

Compound-centric Plot

Use the Compound-centric Plot pane to view an overlay of the quantitation and confirming peaks for each sample in the batch.

Follow these procedures:

- [To select a sample to view](#)
- [To change the target and confirming peaks that are displayed](#)
- [To change the number of rows and columns in the display](#)
- [To copy a plot to another format](#)

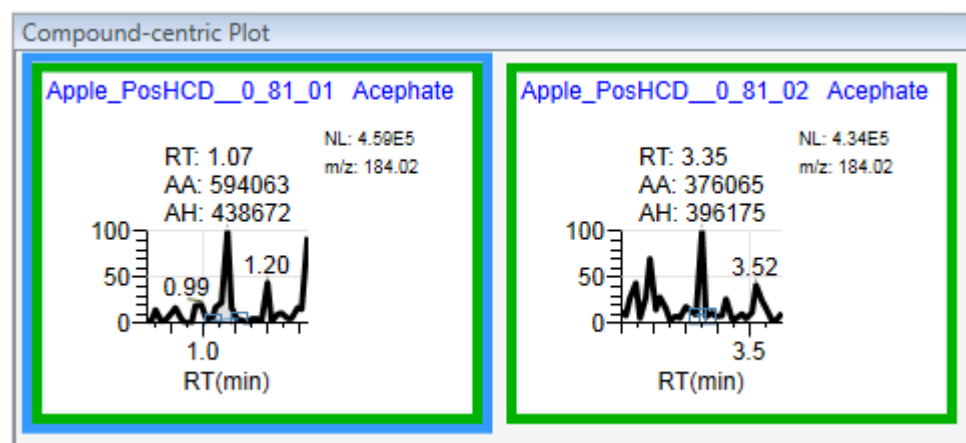
❖ **To select a sample to view**

In the Compounds pane, select the compound that you want to view.

Compounds				Sample Results			
Flags	Compound	RT	Type	Flags	Status	Filename	
1	Acephate	0.00,0.00	Target Compound	1	<input checked="" type="checkbox"/>	2	Apple_PosHCD_0_81
2	Alachlor	0.00,0.00	Target Compound	2	<input checked="" type="checkbox"/>	3	Apple_PosHCD_0_81
3	Ametryn	0.00,0.00	Target Compound	3	<input checked="" type="checkbox"/>	4	Apple_PosHCD_4_05
4	Aminocarb	0.00,0.00	Target Compound				
5	Atrazine	0.00,0.00	Target Compound				

The Compound-centric Plot pane displays the compound chromatograms for the selected compound in each sample in the batch. A blue border indicates the compound in the sample that is currently selected in the Sample Results pane.

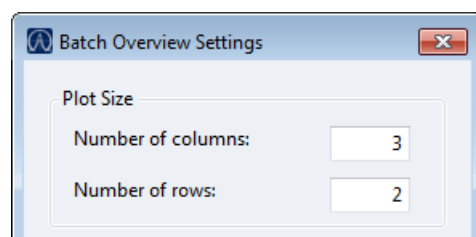
Figure 38. Compound-centric Plot pane



❖ **To change the number of rows and columns in the display**

1. Right-click and choose **Display Settings**.

The Batch Overview Settings dialog box opens.



2. To change the number of rows or columns to fit in the Compound-centric Plot pane, type new values in the Number of Columns or Number of Rows boxes.

4 Using the Analysis Mode for Quantitation Batches

Working in Data Review for Quantitation Batches

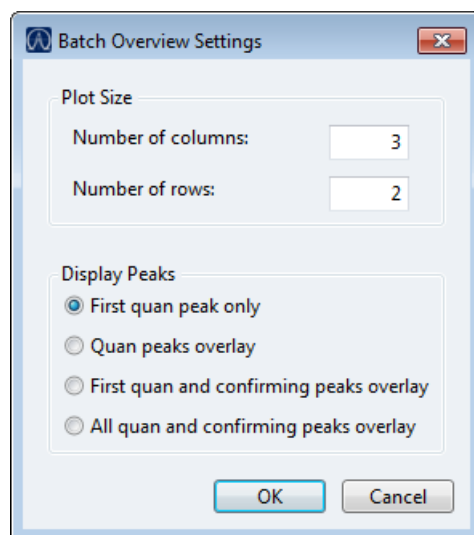
These values do not change the number of rows (compounds) and columns (samples) that are available in the Compound-centric Plot pane. They determine how many rows and columns you want to view at one time in the display. The default is two rows and three columns.

3. Click **OK**.

❖ To change the target and confirming peaks that are displayed

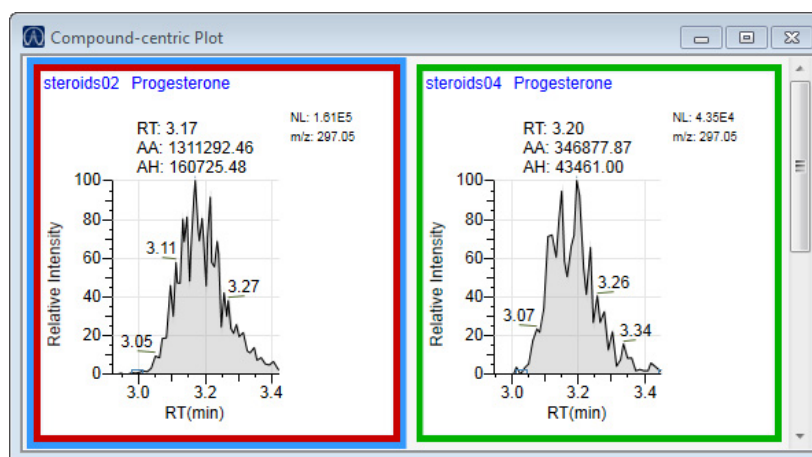
1. Right-click and choose **Display Settings**.

The Batch Overview Settings dialog box opens.

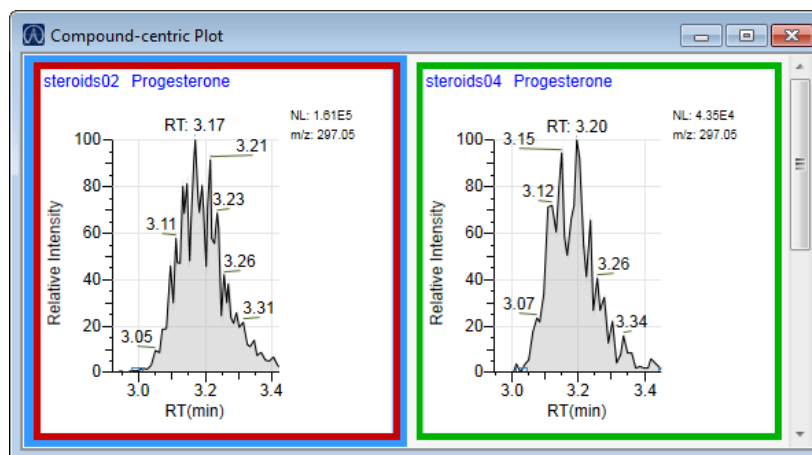


2. Select any of the following display options:

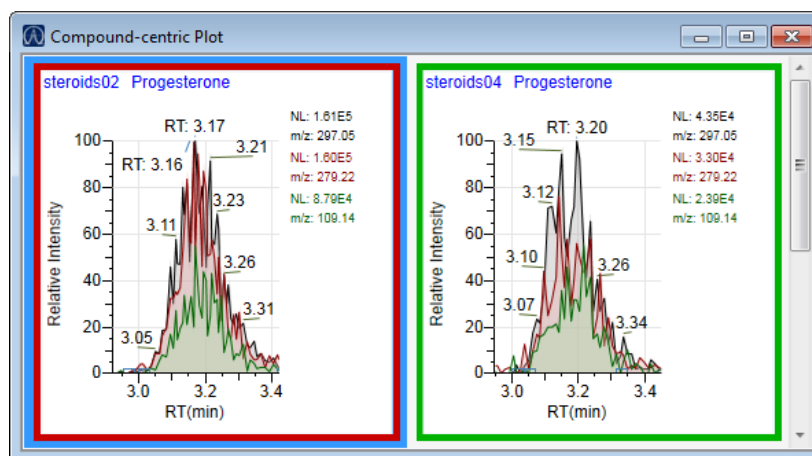
- First Quan Peak Only



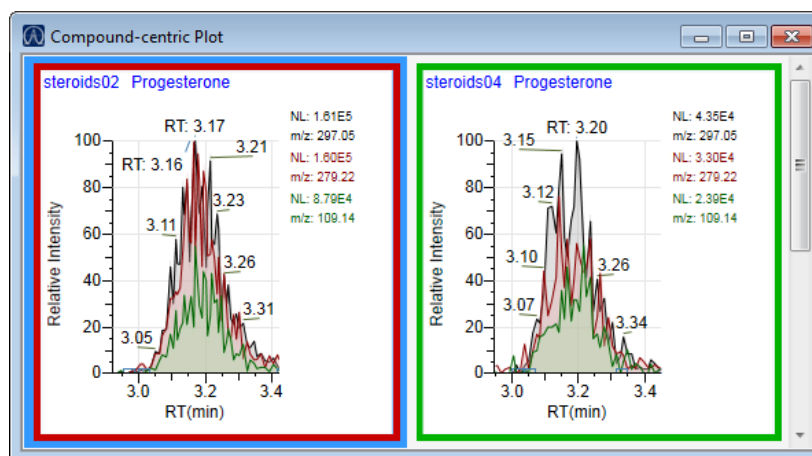
- Quan Peaks Overlay



- First Quan and Confirming Peaks Overlay



- All Quan and Confirming Peaks Overlay



3. Click **OK**.

❖ To zoom in on a peak

1. In any of the chromatogram cells, drag your cursor to delineate a rectangle around the peak.

The delineated area expands to fill the chromatogram area.

2. To restore the default view, right-click the chromatogram plot and choose **Reset Scaling**.

❖ To copy a plot to another format

1. Right-click the plot that you want to copy and choose **Copy to Clipboard**.

The application copies the plot graphic to the Clipboard.

2. In the source format (for example, an Excel spreadsheet or email window), right-click and choose **Paste**.

The application pastes the graphic into the new format.

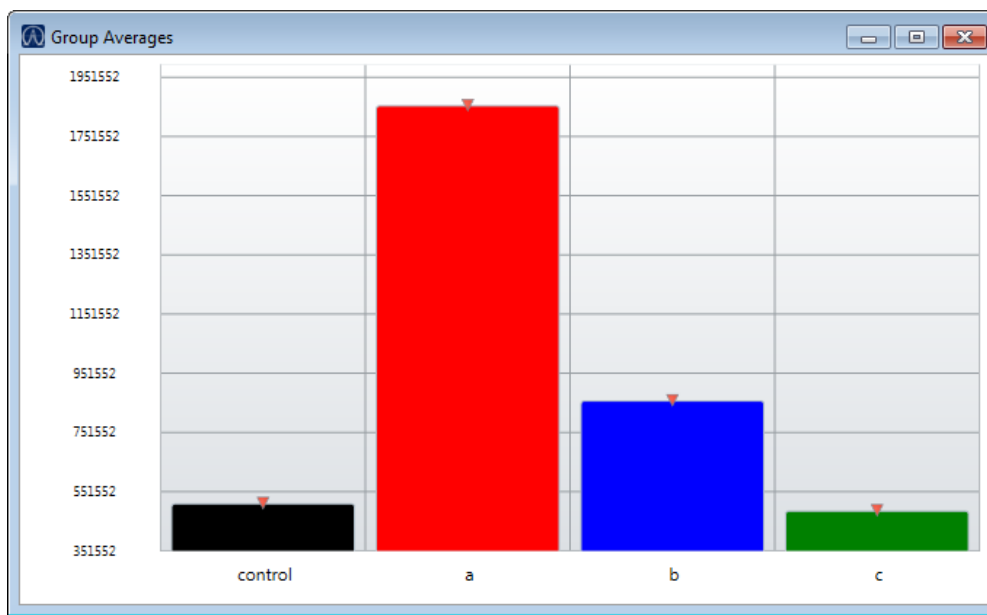
Group Averages

Use the [Group Averages pane](#) to compare the peak areas of different samples to a control group of samples. You must have at least two groups defined, and one group must be a control group.

The control group is displayed to the left, and other samples and groups are listed in the order in which they appear in the Batch View. To specify a group as the control group, see [Groups](#).

Each group displays a colored bar that indicates the average response of the group.

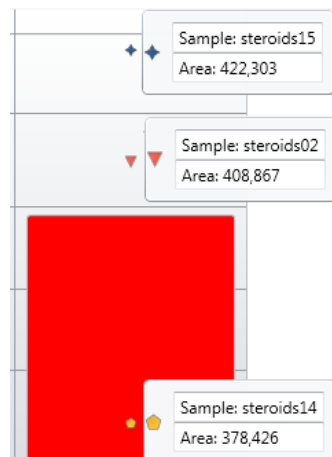
Figure 39. Group Averages pane



❖ To display areas for each sample in the batch

Hold the cursor over an indicator icon (▼, ◆).

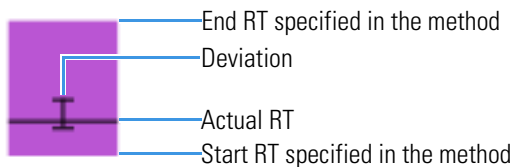
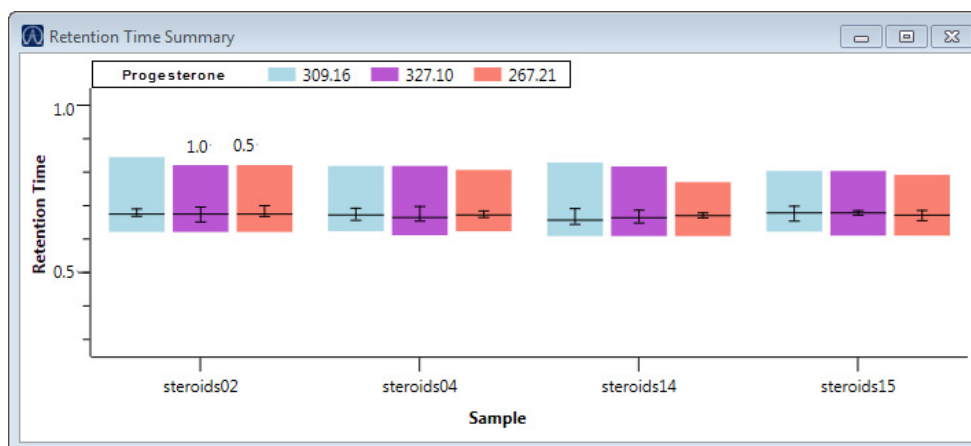
The application displays the area for the selected compound in each represented sample.



Retention Time Summary

Use the Retention Time Summary pane to view variations of the retention times for a compound across all samples in a batch. The Retention Time Summary pane displays each color-coded compound for each sample in the batch.

Figure 40. Retention Time Summary pane



Caution Flags

In the Compound View, the application displays caution flags in both the Compounds pane and in the Sample Results pane.

See the following topics:

- [Flags in the Compounds Pane](#)
- [Flags in the Sample Results Pane](#)
- [Error Indicators in the Compound-centric Plot pane](#)


Flags in the Compounds Pane

The Flags column in the Compounds pane displays a caution flag if the compound in any of the samples is not in compliance with the method criteria.

The screenshot shows two panes: 'Compounds' and 'Sample Results'. In the 'Compounds' pane, the 'Flags' column shows a yellow caution flag for 'Sulfisomidine' (Expected RT: 1.92). In the 'Sample Results' pane, the 'Flags' column shows three rows of data for 'Sulfisomidine' in samples B_25557, B_26154, and B_26473. Each row has a flag icon (green, red, and orange respectively) and a status indicator (green circle). Labels with arrows point to the 'Sulfisomidine' row in the Compounds pane and the flags in the Sample Results pane.

Selected Sulfisomidine compound

Flags for the Sulfisomidine compound in each sample

Click the caution flag icon, , to display the details. Information in the pop-up box shows the sample where the compound is in error and describes the exact error condition.


The screenshot shows the 'Compounds' pane with 'FENTHION-CE20-R20-TL75-QED' selected. A pop-up box is displayed over the 'Flags' column, showing the sample name 'Benzo_25557' and the error condition: '1.909: Apex retention time differential of 0.137 > coelution 0.025'. Labels with arrows point to the sample name and the error condition in the pop-up box.



Sample name

Error condition





Flags in the Sample Results Pane

The Flags column in the Sample Results pane displays a flag if the selected compound in the sample is not in compliance with the method criteria.

Hold your cursor over the flag icon, , to display the details for the selected compound in the sample.

Sample Results					
	Flags	Flag Details	Status	Filename	Sample Type
1		I		B_25557	Specimen
2	1.909: Apex retention time differential of 0.137 > coelution 0.025				

Colored flags in the Sample Results pane have the following meaning.

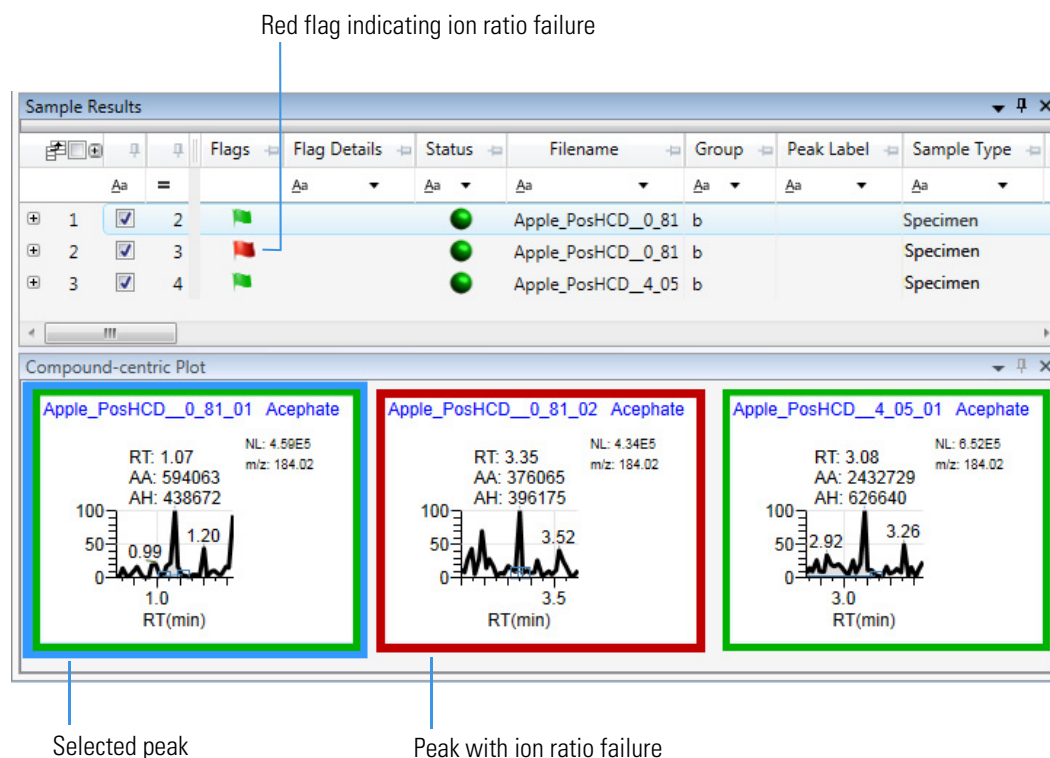
Flag	Description
	Green for compounds that are over the LOR amount specified in the method.
	Red for compounds that have violated (or are activated by) any of the values set in the method. Refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the <i>TraceFinder Lab Director User Guide</i> .
	Red for compounds that are outside the specified ion ratio range. See Ion ratio failure flag .
	Orange for compounds that are not found in Calibrator or QC sample types. “Not found” compounds are below the LOQ, below the LOD, or between the LOD and LOQ values specified in the method. The Sample Results pane does not flag compounds that are not found in Specimen sample types.
	No flag for compounds that have no errors or where no report options are selected.

Note These criteria for flag states do not apply to Negative sample types when the compound is an internal standard.

Error Indicators in the Compound-centric Plot pane

In the Compound-centric Plot pane, peak plots are outlined with the color of their associated error flag. In the following example, the peak plot is highlighted in blue to indicate that Benzo_25557 is the selected sample and outlined in red to indicate that the FENTHION compound in the selected sample is outside the specified ion ratio range.

Figure 41. Ion ratio failure flag



Compound Details for Quantitation Batches

The Compound Details pane is identical for Sample View and Compound View. Use the Compounds Details pane to display any of the following types of data:

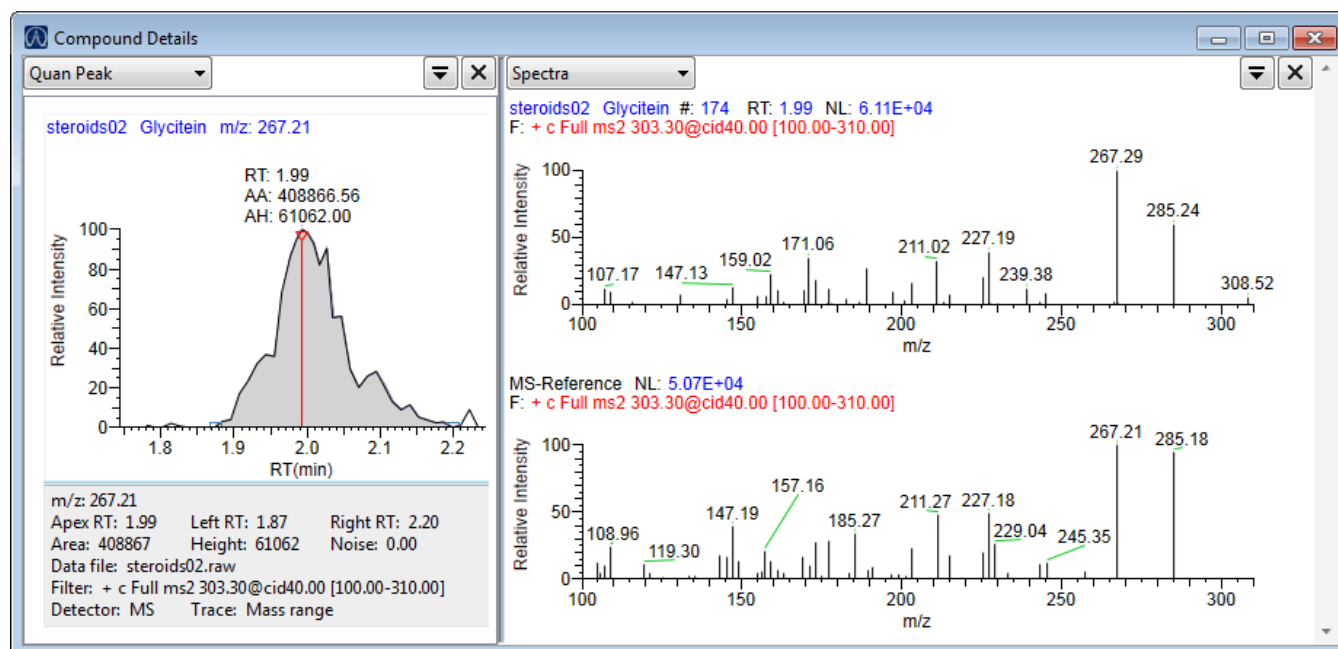
- [Quan Peak](#)
- [Confirming Ions](#)
- [Calibration Curve](#)
- [Ion Overlay](#)
- [ISTD](#)
- [Reference Peak](#)
- [Spectra](#)
- [Library Match](#)

- Isotope
- Fragments
- Quan Peaks Overlay
- Caution Flags

❖ **To open the Compound Details pane**

1. Double-click the chromatogram in the Sample-centric plot or the Compound-centric plot.

The Compound Details pane opens.



By default, the first display pane shows the quantitative peak for the selected compound.

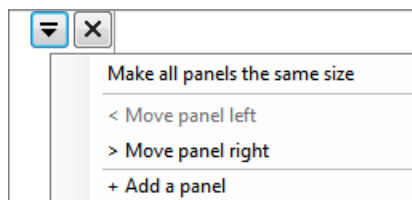
2. In the second pane, select the additional type of data that you want to display.

Follow these procedures to change the display of the peak data in either of the panes:

- To change the peak panes
- To zoom in on a peak

❖ To change the peak panes

In any of the peak panes, click  to view a list of commands.



Command	Description
Make All Panels the Same Size	Evenly divides the area to make all panes the same width. This command does not change the pane height.
Move Panel Left	Moves the current pane one space to the left. This command is not available when the current pane is the leftmost pane.
Move Panel Right	Moves the current pane one space to the right. This command is not available when the current pane is the rightmost pane.
Add a Panel	Adds an empty peak pane to the display. You can display a maximum of four peak panes.

❖ To zoom in on a peak

1. In any of the views, drag your cursor to delineate a rectangle around the peak or spectrum.

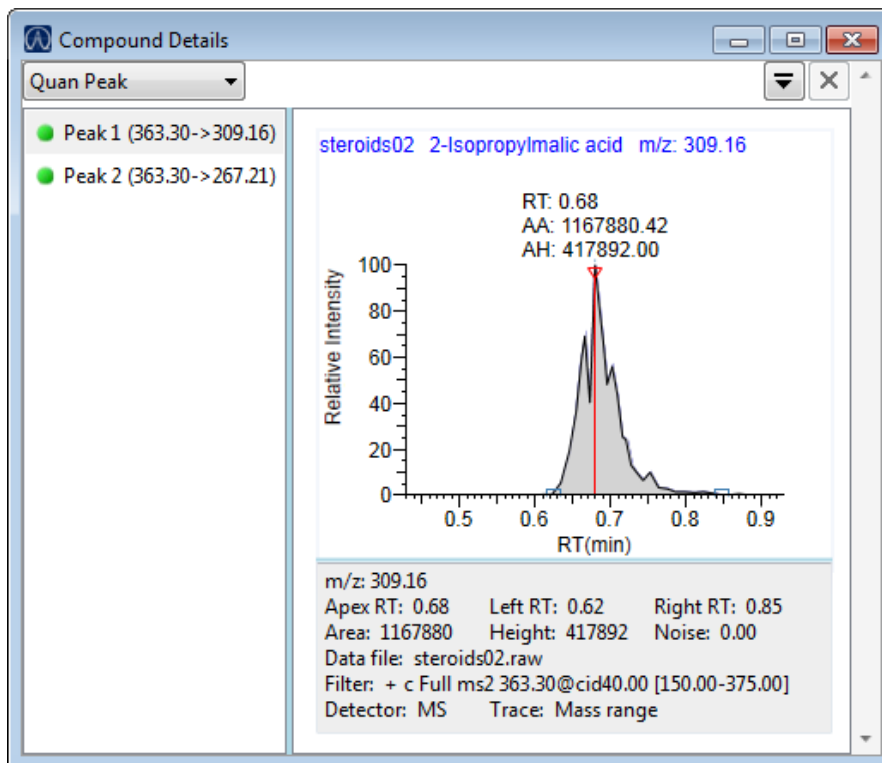
The delineated area expands to fill the view.

2. To restore the default view, right-click the chromatogram plot and choose **Reset Scaling**.

Quan Peak

A compound can have multiple quantitative peaks. You can switch between quantitative peaks, but you cannot view multiple quantitative peaks at the same time. The Quan Peak pane uses a unique shortcut menu. See [Quan Peak pane shortcut menu commands](#).


Figure 42. Quantitative peak pane with multiple quantitative peaks



Follow these procedures to modify the quantitative peak data:

- To manually integrate a peak by swiping the peak
- To manually integrate a peak by moving the peak delimiters
- To remove a manually created peak
- To switch between Method, Manual, and User integration modes
- To customize the labels in the display
- To modify the peak detection settings

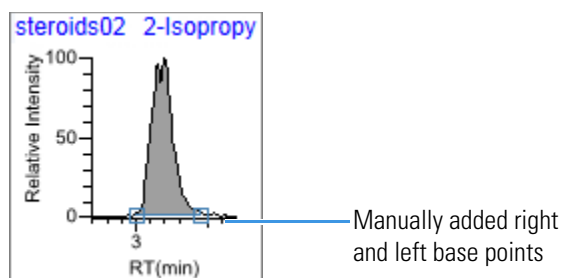
❖ To manually integrate a peak by swiping the peak

1. Right-click anywhere in the quantitative peak pane and then choose **Swipe Manual Integration**, or click the icon in the upper right corner, .
2. Click the left base of the peak that you want to identify.
3. Drag to the right base and release the mouse.

4 Using the Analysis Mode for Quantitation Batches

Working in Data Review for Quantitation Batches

The application places the peak delimiter tags at these locations and automatically updates the peak values (area, height, and so forth) in the result set.



❖ To manually integrate a peak by moving the peak delimiters

1. Hold your cursor over one of the two peak delimiter tags in the peak pane.

When the tag can be selected, the cursor changes to a crosshair-style cursor. You can zoom in on the baseline to make it easier to select the tag.



2. Drag the peak delimiter tag to another location and automatically update the peak values (area, height, and so forth) into the result set.

The generated reports for the data identify the manual modifications.

You can store two peak value sets (method and manual integration settings) with each compound in each file. These settings can result in a different set of stored values. The application originally calculates the method values based on the processing method parameters. The manual values are a result of what you edited.

❖ To remove a manually created peak

Right-click the pane and then choose **Delete Peak**, or click the icon in the upper right corner, .

The application removes the manual integration from the selected peak.

❖ To switch between Method, Manual, and User integration modes

Right-click the chromatogram pane and choose **Integration Settings**.

Initially, the method, manual, and user integration settings that are stored for a compound and file are identical. When you switch modes, the saved result set does not change. However, when either manual or user data are available, both the chromatogram plots and the result set update as you switch between modes.

As you switch between modes, each pane in Data Review reflects the changes. The generated reports for the data identify the manual or user modifications.

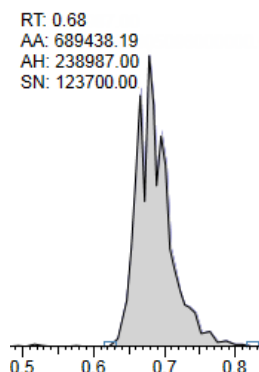
❖ **To customize the labels in the display**

1. Right-click and choose **Peak Display Settings**.

The **Peak Display Settings** dialog box opens.

2. To display labels for the peak retention time (RT), area (AA), height (AH), or signal-to-noise (SN), select the appropriate check box.
3. Click **OK** in the Peak Display Settings dialog box.

Figure 43. Example of displayed labels

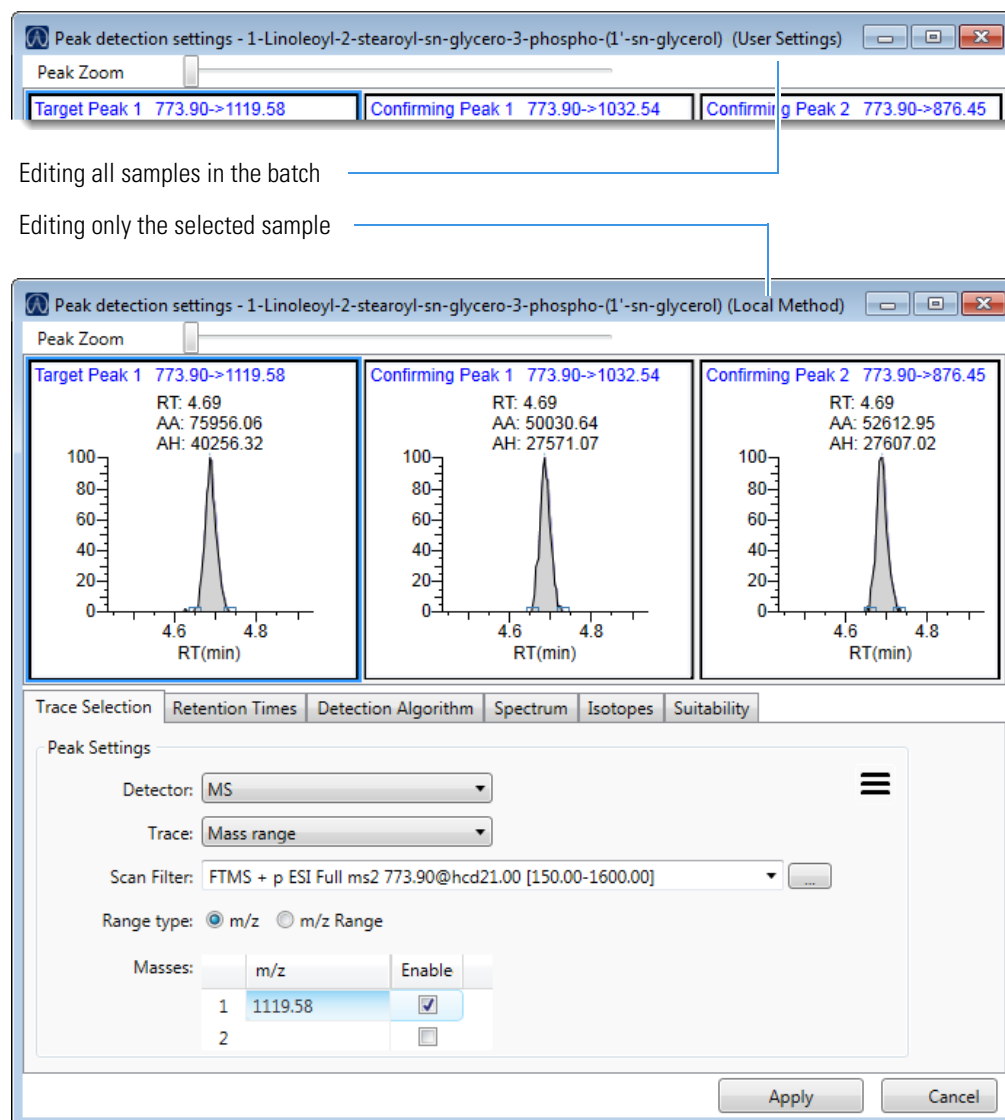


❖ **To modify the peak detection settings**

1. Right-click the chromatogram view and choose one of the following:
 - **Peak Detection Settings > Edit Local Method Peak Detection Settings:** Makes changes to the selected compound for all samples in this batch. The application identifies these settings as “Manual.”
 - **Peak Detection Settings > Edit User Peak Detection Settings:** Makes changes to the selected compound for only the selected sample. The application saves these changes with the batch and stops applying the local method detection settings to the compound for this sample only. The application identifies these settings as “User.”

The Peak Detection Settings dialog box opens where you can adjust detection settings that were specified in the method. The title bar of the dialog box lists the selected compound and indicates whether you are making changes to only the selected sample or to the local method.

Figure 44. Peak Detection Settings dialog box





2. Edit any of the detection settings.

For detailed descriptions of all detection settings, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

3. To save your changes to this compound, click **Apply**.
 - When you are editing a single sample, the application makes changes to the selected compound for this sample. If the sample is a calibration sample type, this update changes the calibration curve which, in turn, affects all calculated amounts.
 - When you are editing the local method, the application makes changes to the selected compound for all samples in this batch.

Table 30. Quan Peak pane shortcut menu commands

Command	Description
Reset Scaling	Resets the original scaling after a zoom operation.
Copy to Clipboard	Copies the graphic display to the Clipboard.
Send RT to Method	Sets the current retention time as the expected retention time for the compound in the local method.
Swipe Manual Integration	Switches to manual integration mode, where you can manually define the peak. Also available from the icon in the upper right corner,  .
Integration Settings (Method, Manual, User)	Switches the display between Method, Manual, or User integration modes. Method (default): The settings defined in the master method. Manual: Modifications made using the Peak Detection Settings > Edit Local Method Peak Detection Settings command. The application applies the modified peak detection settings to all samples in the batch. User: Modifications made using the Peak Detection Settings > Edit User Peak Detection Settings command. The application applies the modified peak detection settings to only the selected sample.
Clear User Peak Detection Settings	Resets modified peak detection settings to the settings in the master method, and identifies the integration mode as Method.
Delete Peak	Removes the peak and all data for this peak from all Compound View panes. Also available from the icon in the upper right corner,  .
Peak Display Settings	Opens the Peak Display Settings dialog box .
Peak Detection Settings	Edit Local Method Peak Detection Settings: Edits the peak detection settings for all samples in the batch. The application identifies these settings as “Manual.” Edit User Peak Detection Settings: Edits the peak detection settings for only the selected sample. The application identifies these settings as “User.” After you apply either of these updates, the application does not retain the modified integration settings. The changes that you make on the Signal page in the Peak Detection Settings dialog box can affect all quantification and ion ratio calculations and are not auto-reversible. These changes can also create a mismatch with calibration files that are not fully defined in the batch.
Note The Peak Detection Settings commands are also available in the Confirming Ions pane.	

Confirming Ions

The Confirming Ions pane displays a list of peaks, a graphical view of all qualifying/confirming ions for the selected compound, and calculated ion ratios and ion ratio acceptance windows. A red border indicates that an ion ratio is outside of its window. See [Quantitative peak with multiple confirming ions](#). The Confirming Ions pane uses a unique shortcut menu. See [Confirming Ions pane shortcut menu commands](#).

In the peaks list, the application displays the mass of each confirming ion peak. When a peak is within the tolerance set, the peak status is green. When a peak is not within the tolerance set, the peak status is red.

Note For compounds with an analog detection type, the application displays “No Confirming Ions are Enabled” in the Confirming Ions pane.

Follow these procedures:

- [To remove a peak](#)
- [To customize the labels in the display](#)
- [To display surrounding apex times](#)
- [To manually integrate a peak](#)
- [To switch between Method, Manual, and User integration modes](#)

❖ To remove a peak

Right-click the Confirming Ions pane, and choose **Delete Peak**.

The application removes the peak displayed in the Confirming Ions pane. All data for this peak are removed from the Compound View panes.

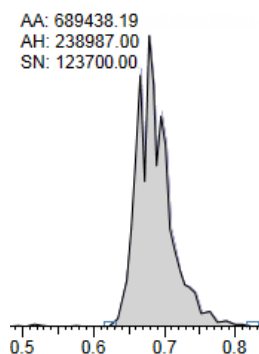
❖ To customize the labels in the display

1. Right-click and choose **Peak Display Settings**.

The [Peak Display Settings Dialog Box](#) opens.

2. To display labels for the peak retention time (RT), mass area (MA), mass height (MH), or signal-to-noise (SN), select the appropriate check box. See [Example of displayed labels](#).
3. Click **OK** in the Peak Display Settings dialog box.

Figure 45. Example of displayed labels



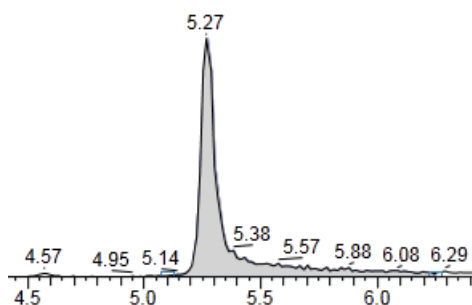
❖ **To display surrounding apex times**

1. Right-click and choose **Peak Display Settings**.

The **Peak Display Settings** dialog box opens.

2. Select the **Show All Apex Times** check box.
3. Click **OK** in the Peak Settings dialog box.

Figure 46. Example of displayed apex times



❖ **To manually integrate a peak**

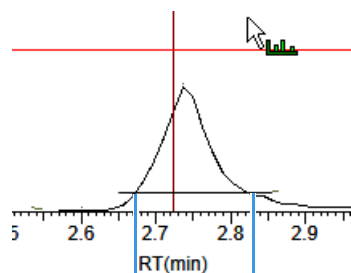
1. Right-click and then choose **Swipe Manual Integration**, or click the **Swipe Manual Integration** icon, .

The cursor changes to indicate the manual integration mode.

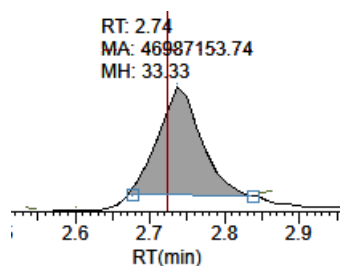


2. Drag the cursor to define the beginning and ending base points for the new peak.

Note You must drag the cursor inside the x axis and y axis.



Drag the cursor across the peak at the base points for the new peak.



The application identifies the peak and indicates the manual integration in the labels.

The application updates the area of the peak to the baseline that you have just added. When you regenerate the reports, they display the manually modified data.

You can store three peak value sets (method, manual, and user integration settings) with each compound in each file. These settings can result in a different set of stored values. The application originally calculates the method values based on the processing method parameters. The manual or user values are a result of what you edited.

Note Because a Blank Report displays only the quantitation mass, when you manually integrate a confirming ion, the manual integration flag in the report is displayed on the quantitation mass.

❖ To switch between Method, Manual, and User integration modes

Right-click the Confirming Ions pane and choose **Integration Settings**.

Initially, the method, manual, and user integration settings that are stored for a compound and file are identical. When you switch modes, the saved result set does not change. However, when either manual or user data are available, both the chromatogram plots and the result set update as you switch between modes.

As you switch between modes, each pane in Data Review reflects the changes. The generated reports for the data identify the manual or user modifications.

Figure 47. Quantitative peak with multiple confirming ions

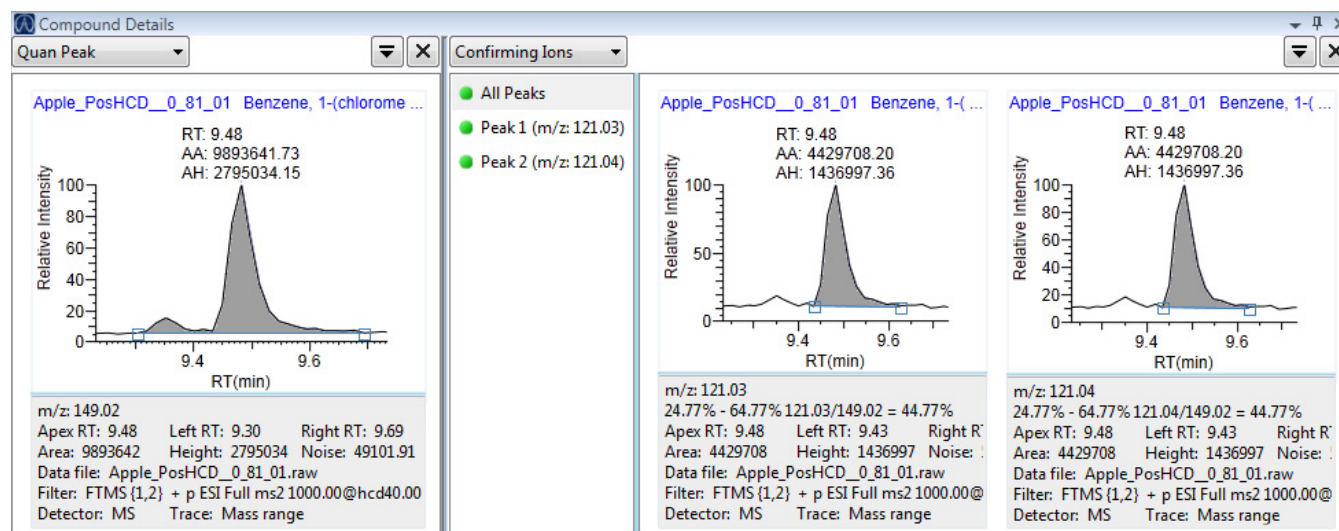


Figure 48. Confirming ion with coelution failure

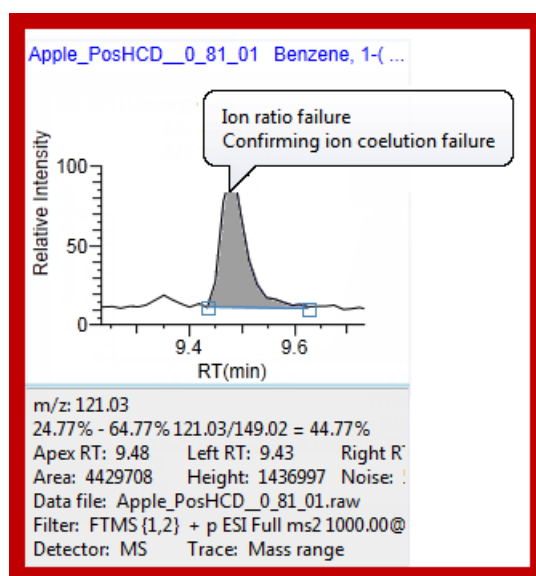


Table 31. Confirming Ions pane shortcut menu commands (Sheet 1 of 2)



Command	Description
Reset Scaling	Resets the original scaling after a zoom operation.
Copy to Clipboard	Copies the graphic display to the Clipboard.
Swipe Manual Integration	Switches to manual integration mode, where you can manually define the peak.
Also available from the icon in the upper right corner,  .	

Table 31. Confirming Ions pane shortcut menu commands (Sheet 2 of 2)

Command	Description
Integration Settings (Method, Manual, User)	<p>Switches the display between Method, Manual, or User integration modes.</p> <p>Method (default): The settings defined in the master method.</p> <p>Manual: Modifications made using the Peak Detection Settings > Edit Local Method Peak Detection Settings command. The application applies the modified peak detection settings to all samples in the batch.</p> <p>User: Modifications made using the Peak Detection Settings > Edit User Peak Detection Settings command. The application applies the modified peak detection settings to only the selected sample.</p>
Clear User Peak Detection Settings	Resets modified peak detection settings to the settings in the master method, and identifies the integration mode as Method.
Delete Peak	<p>Removes the peak and all data for this peak from all Compound View panes.</p> <p>Also available from the icon in the upper right corner, .</p>
Peak Display Settings	Opens the Peak Display Settings dialog box .
Peak Detection Settings	<p>Edit Local Method Peak Detection Settings: Edits the peak detection settings for all samples in the batch. The application identifies these settings as “Manual.”</p> <p>Edit User Peak Detection Settings: Edits the peak detection settings for only the selected sample. The application identifies these settings as “User.”</p> <p>After you apply either of these updates, the application does not retain the modified integration settings.</p> <p>The changes that you make on the Signal page in the Peak Detection Settings dialog box can affect all quantification and ion ratio calculations and are not auto-reversible. These changes can also create a mismatch with calibration files that are not fully defined in the batch.</p>
<p>Note The Peak Detection Settings commands are also available in the Confirming Ions pane.</p>	

Peak Display Settings Dialog Box

Use the features in the Peak Display Setting dialog box to change the labels in the Confirming Ions pane display.

Figure 49. Peak Display Settings dialog box

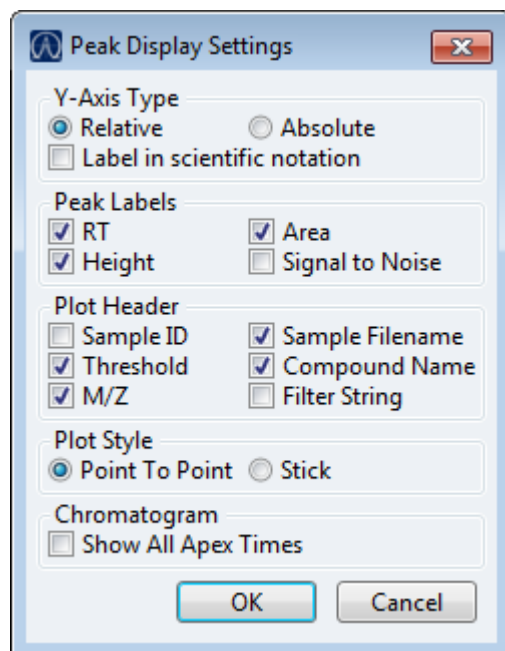
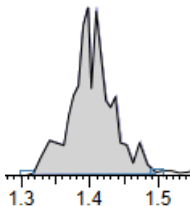
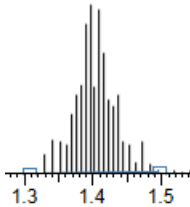


Table 32. Peak Display Settings dialog box parameters (Sheet 1 of 2)

Command	Description
Y-Axis Type	
Relative	Displays the <i>y</i> -axis scale from 0 through 100.
Absolute	Displays the <i>y</i> -axis scale from 0 to the actual value of the most intense peak in the group.
Label in Scientific Notation	Displays the <i>y</i> -axis scale in scientific notation.
Peak Labels	
RT	Displays the retention time (RT) label in the peak chromatogram pane.
Height	Displays the peak height (AH) label in the peak chromatogram pane.
Area	Displays the peak area (AA) label in the peak chromatogram pane.
Signal to Noise	Displays the signal-to-noise (SN) label in the peak chromatogram pane.
Plot Header	
Sample ID	Displays a user-defined, alphanumeric string that identifies a sample.
Threshold	Displays the peak threshold.

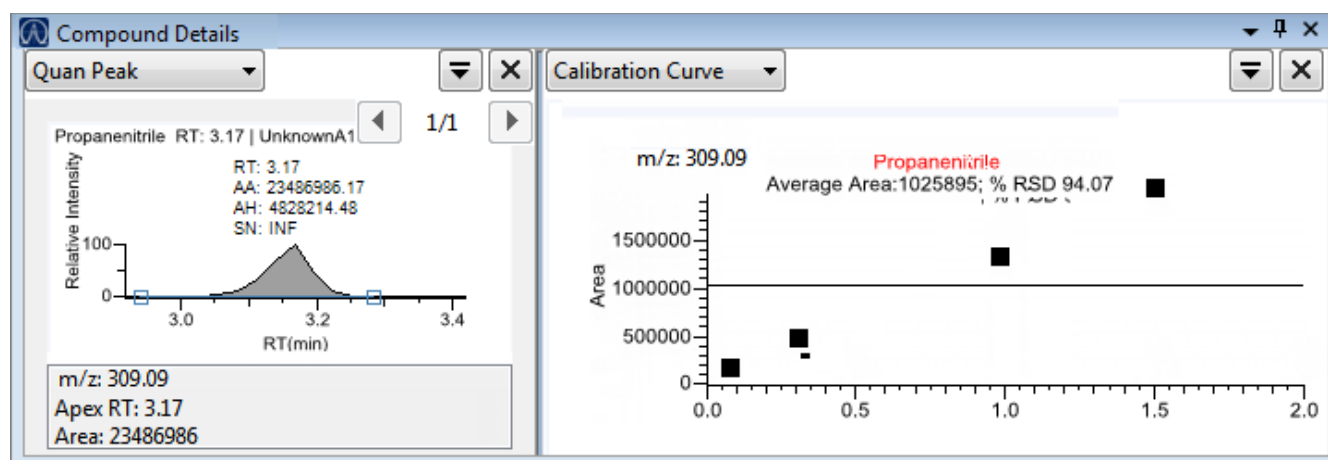
Table 32. Peak Display Settings dialog box parameters (Sheet 2 of 2)

Command	Description
<i>m/z</i>	Displays the mass-to-charge ratio found in the spectrum for the peak. Assumes the charge is 1.
Sample Filename	Displays the name of the current, selected sample where the peak was found.
Compound Name	Displays the matched compound name found in the library or compound database.
Filter String	Displays the filter used to identify the peak. Specified in the raw data file or the master method.
Plot Style	
Point To Point	Displays chromatograms in the following format:
	
Stick	Displays chromatograms in the following format:
	
Chromatogram	
Show All Apex Times	Displays all surrounding apex times in the plot.

Calibration Curve

The Calibration Curve pane displays a graphical view of the calibration curve for the selected compound and key statistical values for evaluating the quality of the calibration. The Calibration Curve pane uses a unique shortcut menu. See [Calibration Curve pane shortcut menu commands](#).

Figure 50. Quantitative peak with a calibration curve plot



❖ To manually exclude a calibration point

In the Sample Results pane, select the **Excluded** check box for the sample.

You might need to use the horizontal scroll bar at the bottom of the table to scroll to the Excluded column. See [Inactive and Excluded Compounds](#).

When the application no longer uses a calibration point, its value appears as an empty box in the graphical view of the calibration curve.

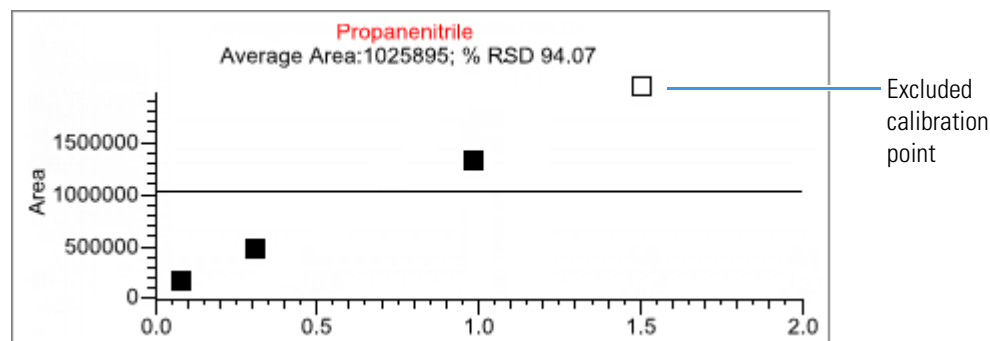


Table 33. Calibration Curve pane shortcut menu commands

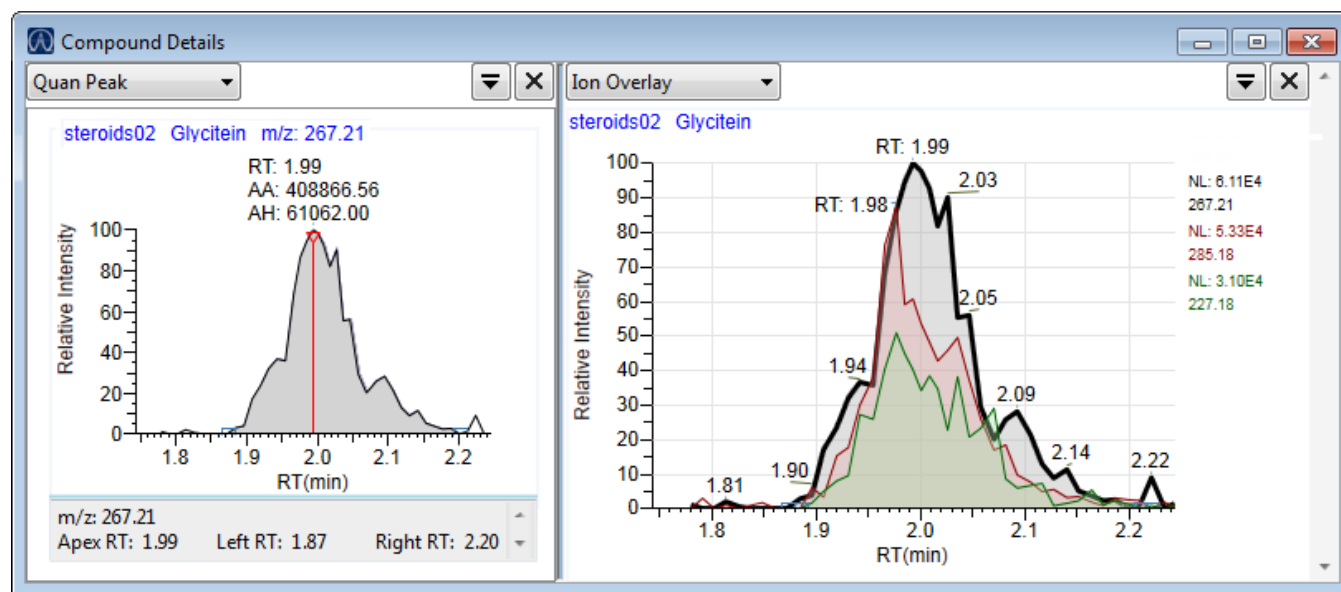
Command	Description
Standard Type	Sets the standard type to External or Internal.
Calibration Curve Type	Sets the calibration curve type to one of the following: <ul style="list-style-type: none"> • Linear: Uses all settings with this exception: When Origin is set to Include, all Weighting values are grayed out and Weighting is set to Equal. • Quadratic: Uses all settings with this exception: When Origin is set to Include, all Weighting values are grayed out and Weighting is set to Equal. • Average RF: Uses no Weighting or Origin selections. All Weighting and Origin values are grayed out. Weighting is set to Equal, and Origin is set to Ignore.
Response Via	Sets the response to Area or to Height.
Weighting	Sets the weighting to equal, 1/X, 1/X ² , 1/Y, or 1/Y ² .
Origin	Sets the origin to Ignore, Force, or Include.
Units	Sets the units.
Done with Settings	Saves the calibration curve settings.
Reset Scaling	Resets the original scale in the calibration curve pane.

Ion Overlay

The Ion Overlay pane represents an overlay of the entire ion set—target and confirming—for the selected compound. Use this pane to graphically review the peak apex alignment and coeluting peak profiles.

Note For compounds with an analog detection type, the application displays “No Data” in the Ion Overlay pane.

Figure 51. Target peak with a confirming ion overlay



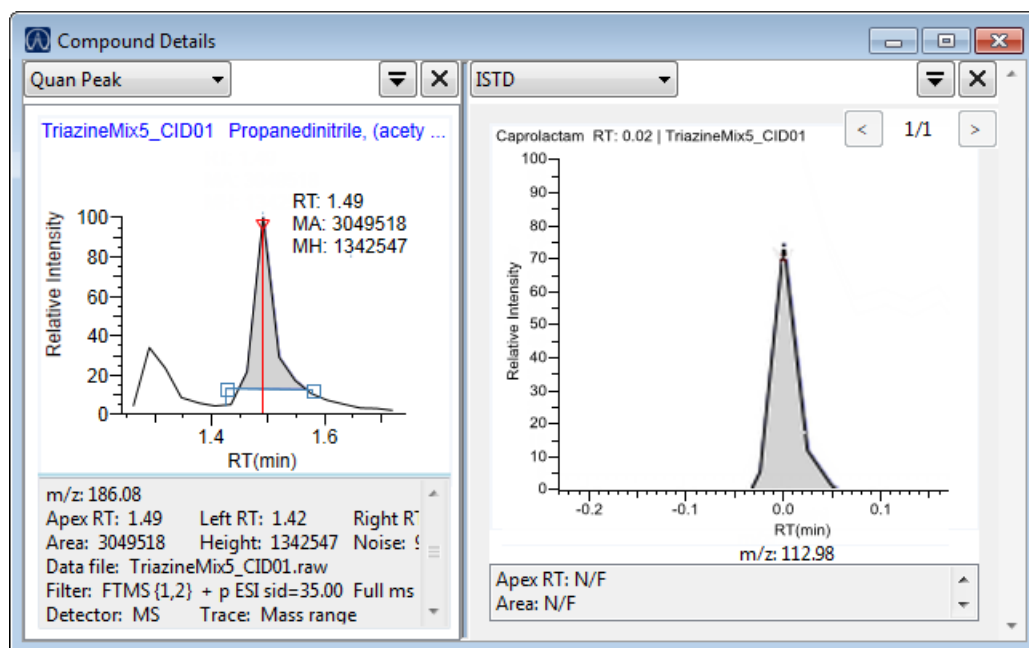
4 Using the Analysis Mode for Quantitation Batches

Working in Data Review for Quantitation Batches

ISTD

The ISTD pane displays the internal standard specified for the compound in the method. Refer to the instructions “To specify an internal standard type for a compound” in Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

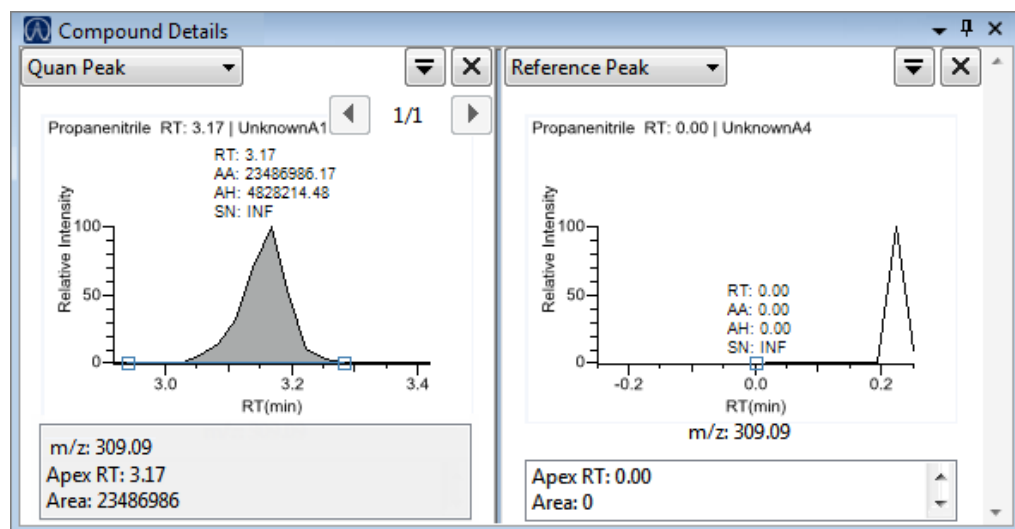
Figure 52. Quantitative peak with an internal standard



Reference Peak

The Reference Peak pane displays the reference peak as specified in the method.

Figure 53. Quantitative peak with a reference peak



❖ **To specify a chromatogram reference peak**

1. In the Batch View task pane, click **Reference Sample**.

An empty reference sample table opens.

Batch View - Batch_A [Quan]


Local Method: Method_A_1

Update

Status	Filename	Sample ID	Sample name	Comment	Vial position	Barcode Actual
--------	----------	-----------	-------------	---------	---------------	----------------

2. Right-click and choose **Add Reference Sample**.

The Open Chromatogram Reference Sample dialog box opens.



Open Chromatogram Reference Sample

C:\TraceFinderData\...\Projects

Project A

Batch_3

UnknownA2

UnknownA3

UnknownA4

UnknownA1

Select a reference sample from the right panel and click "Open".

Open Cancel

Note If you are creating a new method, you will not see any reference samples here. You must create and save a batch using the current method to see the reference samples in this list.

3. Select a project from the list of projects.
4. Select a subproject from the list of subprojects.
5. Select a batch from the list of batches.

The application displays only batches that were created using the current master method.

6. Select a sample from the list of processed samples.

The application displays all the processed samples in the selected batch. To use a sample as a reference sample, the sample must have been processed with the current master method.

7. Click **Open**.

The selected sample is displayed as the chromatogram reference sample in the Method View in the Method Development mode.

Tip To clear the reference sample from the master method, right-click and choose **Delete Selected**.

Spectra

The Spectra pane displays a comparison of the spectra found in the data and the method reference.

Note For compounds with an analog detection type, the application displays “Not Available” in the Spectra pane.

Follow these procedures:

- [To zoom in on a peak](#)
- [To copy a plot to another format](#)
- [To display the MS reference](#)
- [To customize the labels in a small molecule or peptide spectrum](#)

❖ To zoom in on a peak

1. In the Spectra pane, drag the cursor to delineate a rectangle around the peak.

The delineated area expands to fill the view.

2. To restore the default view, right-click the Spectra pane and choose **Reset Scaling**.

❖ To copy a plot to another format

1. In the data spectrum or the reference spectrum, right-click and choose **Copy to Clipboard**.

The application copies the specified plot graphic to the Clipboard.

Note The application copies only the indicated data spectrum or the reference spectrum. It does not copy the entire Spectra pane.

2. In the source format (for example, an Excel spreadsheet or email window), right-click and choose **Paste**.

The application pastes the graphic into the new format.

❖ To display the MS reference

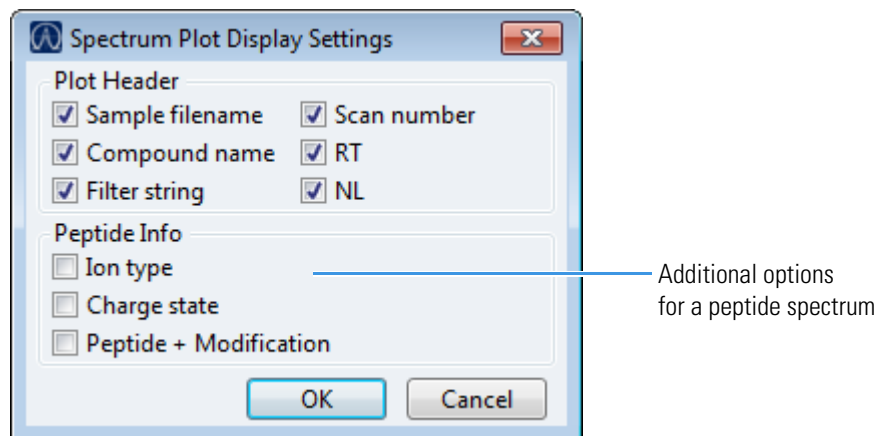
Right-click the Spectra pane and choose **Show MS-Reference**.

The application displays or hides the reference spectrum in the Spectra pane.

❖ To customize the labels in a small molecule or peptide spectrum

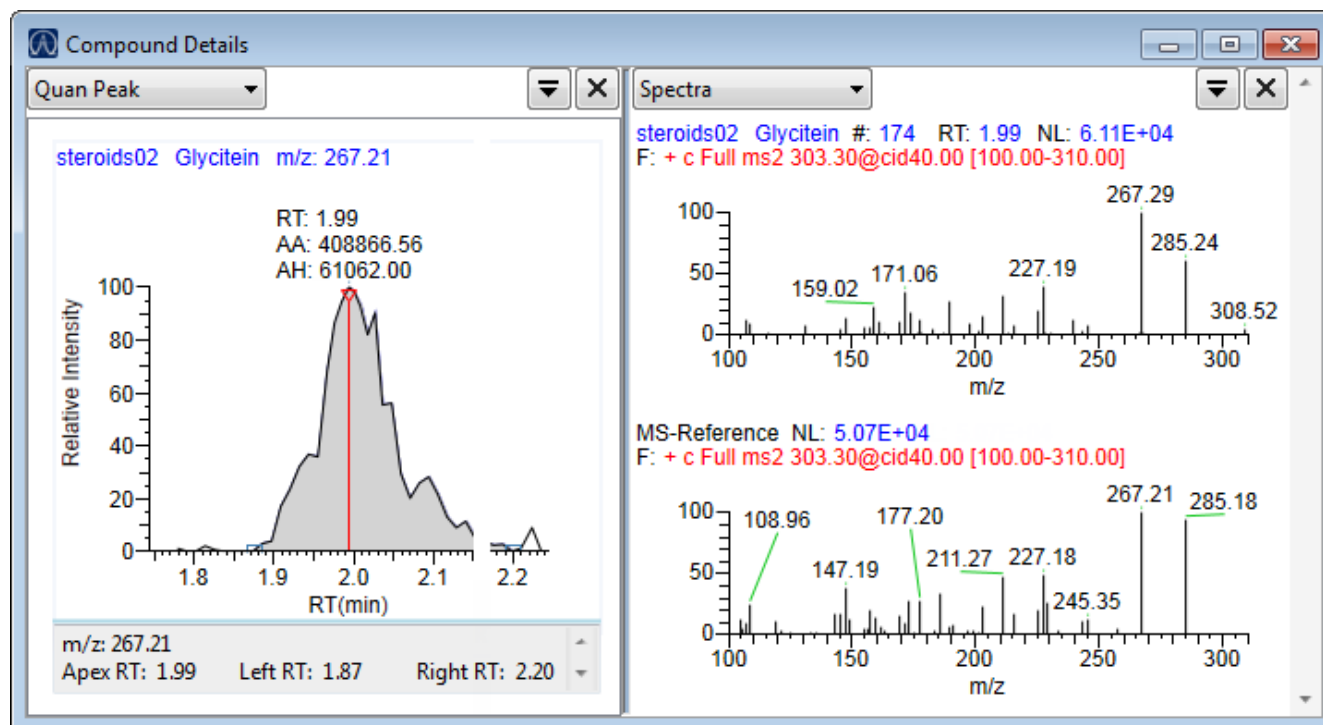
1. Right-click the Spectra pane and choose **Display Settings**.

The Spectrum Plot Display Settings dialog box opens.



2. Select the check box for each label that you want to display and click **OK**.

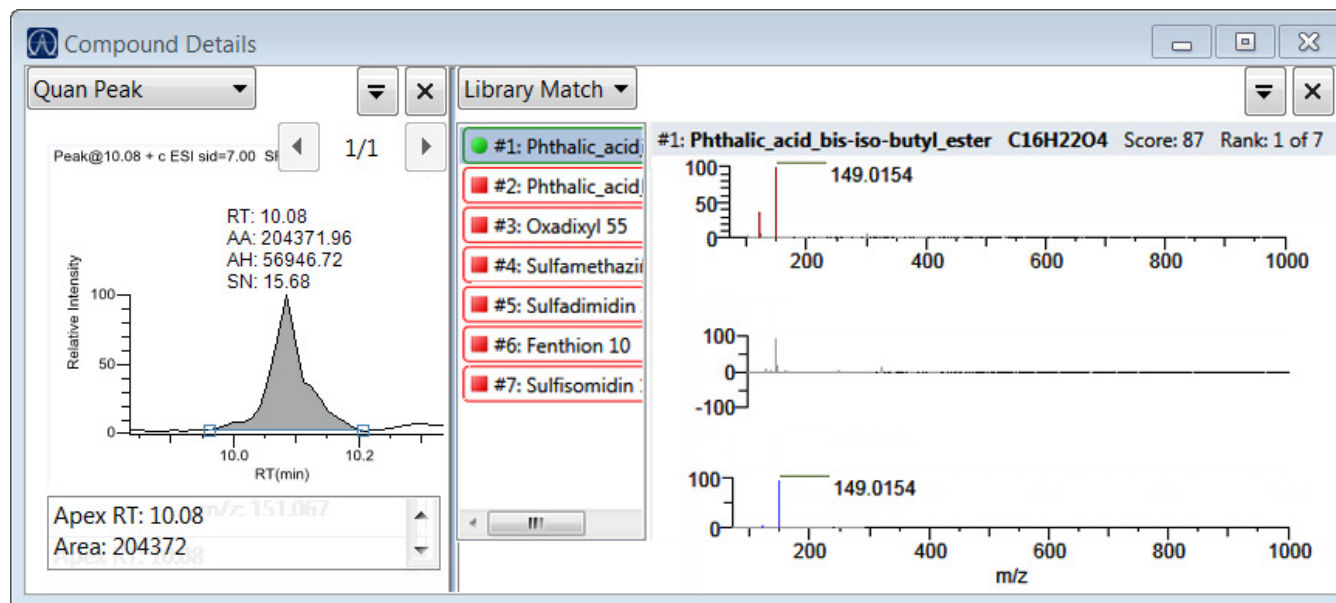
Figure 54. Quantitative peak with data and reference spectra



Library Match

The Library Match pane displays all library matches for the selected compound.

Figure 55. Library Match for the selected compound



If you have no matches for any of your compounds, make sure you have completed all of the following:

- Installed a library.
- Selected screening libraries in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.
- Enabled Library Matching in the method. Refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

Isotope

The [Isotope pane](#) displays all isotopes for the selected compound. The Isotope pane includes a shortcut menu. See [Isotope pane shortcut menu commands](#).

Note When you select different samples, the current compound remains selected (as long as that compound is found in the sample).

The isotopes pane displays isotopic pattern results for all adducts of a compound according to the threshold and deviation parameters defined in the method.

To identify or confirm the presence of a compound, the resulting score percentage from isotopic pattern matching must be higher than the specified fit threshold percentage.

- An isotope peak is not found if its intensity, relative to the monoisotopic ion's intensity, is more than the specified intensity deviation percentage away from the theoretical relative intensity of the isotope ion.
- An isotope peak is found if its measured m/z is less than the specified mass deviation amount away from its expected m/z .

To specify isotopic criteria in a method, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

Figure 56. Isotope pane

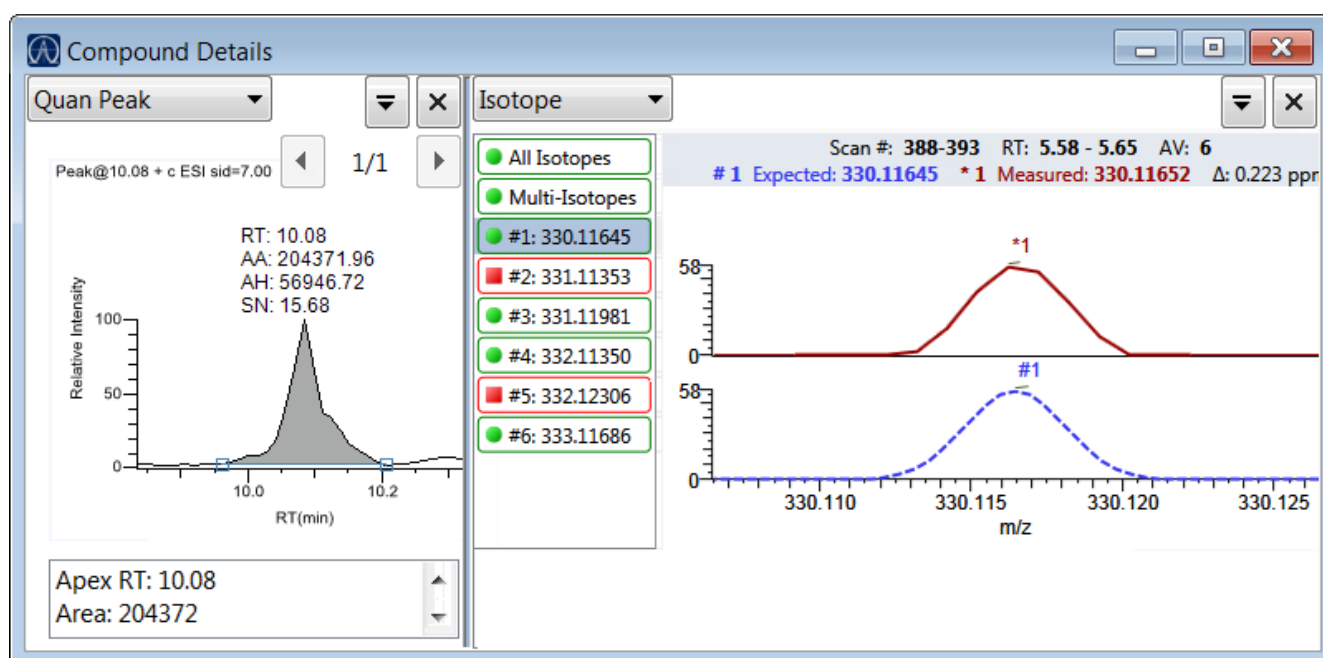


Table 34. Isotope pane shortcut menu commands

Command	Description
Reset Scaling	Resets the original scaling after a zoom operation.
Copy to Clipboard	Copies the graphic display to the Clipboard.
Display Overlay Spectra	Overlays the two spectrum displays, or stacks the peak apex spectrum over the simulated spectrum.
Display Stack Spectra	Overlays the two spectrum displays, or stacks the peak apex spectrum over the simulated spectrum.
Show/Hide Noise Label	Adds a noise label to each peak. Expected isotope peaks (displayed in blue) do not display a noise label.
Show/Hide Resolution Label	Adds a resolution label to each peak. Expected isotope peaks (displayed in blue) do not display a resolution label.

Fragments

The Fragments pane displays fragments in one of two ways:

- All Fragments
- Individual Fragments

All Fragments

The All Fragments view displays a composite of all fragments found in the peak. The application displays the measured peak as a red line; the application displays the expected peak as a blue line.

The application displays these headers for the All Fragments view:

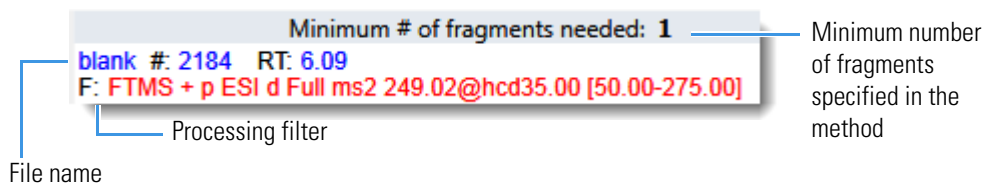


Figure 57. Fragments pane with overlaid spectra for all fragments

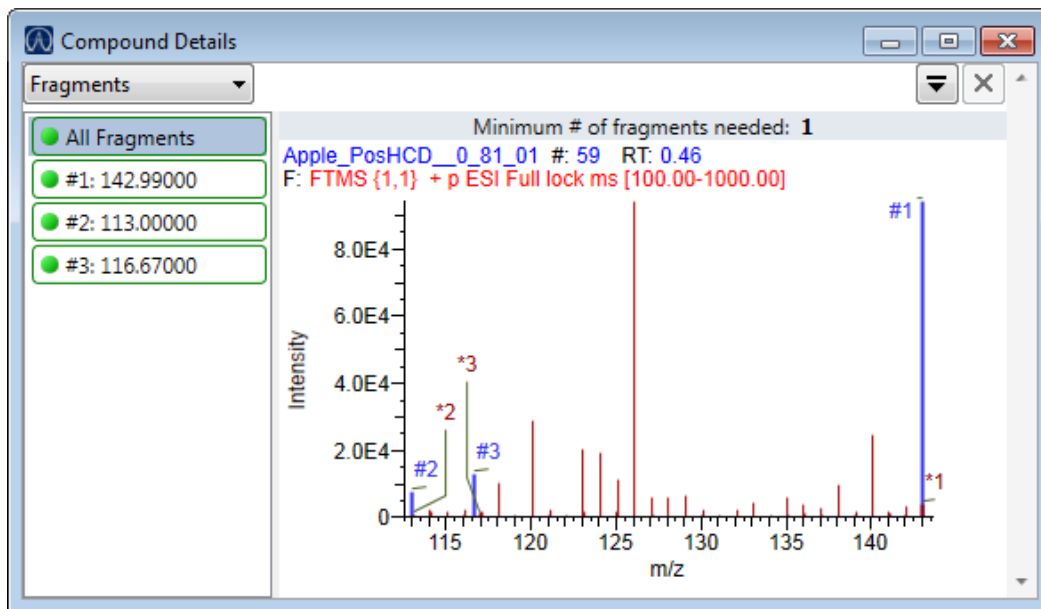
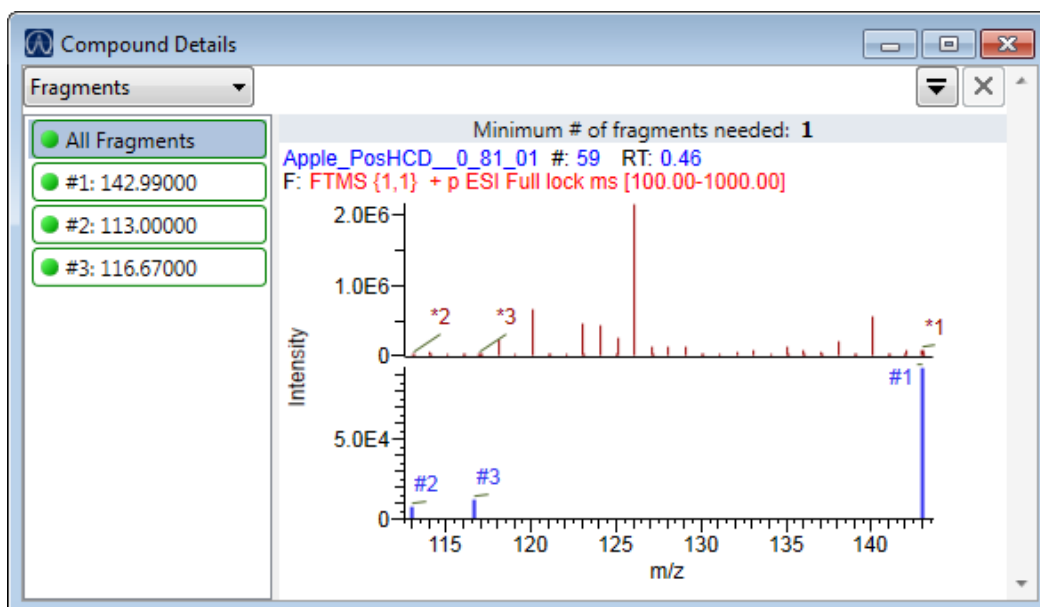


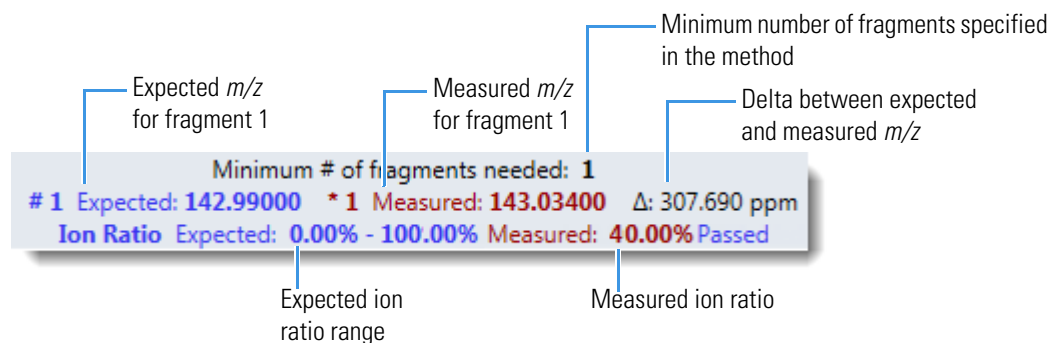
Figure 58. Fragments pane with stacked spectra for all fragments



Individual Fragments

The individual fragments view displays the expected and measured peaks for a single fragment.

The application displays these headers for the individual fragments view:



4 Using the Analysis Mode for Quantitation Batches

Working in Data Review for Quantitation Batches

Figure 59. Fragments pane with overlaid spectra for a single fragment

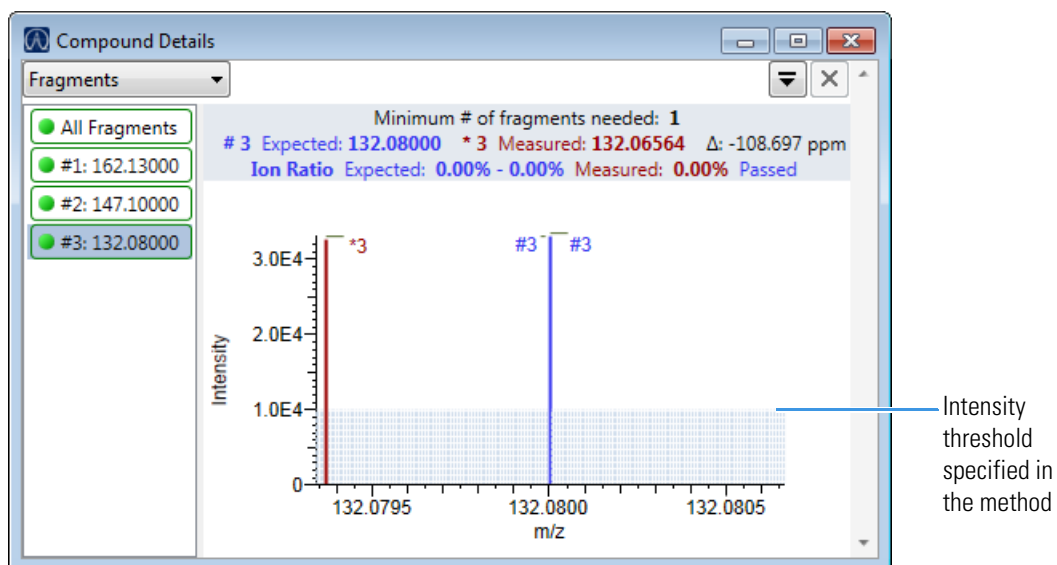
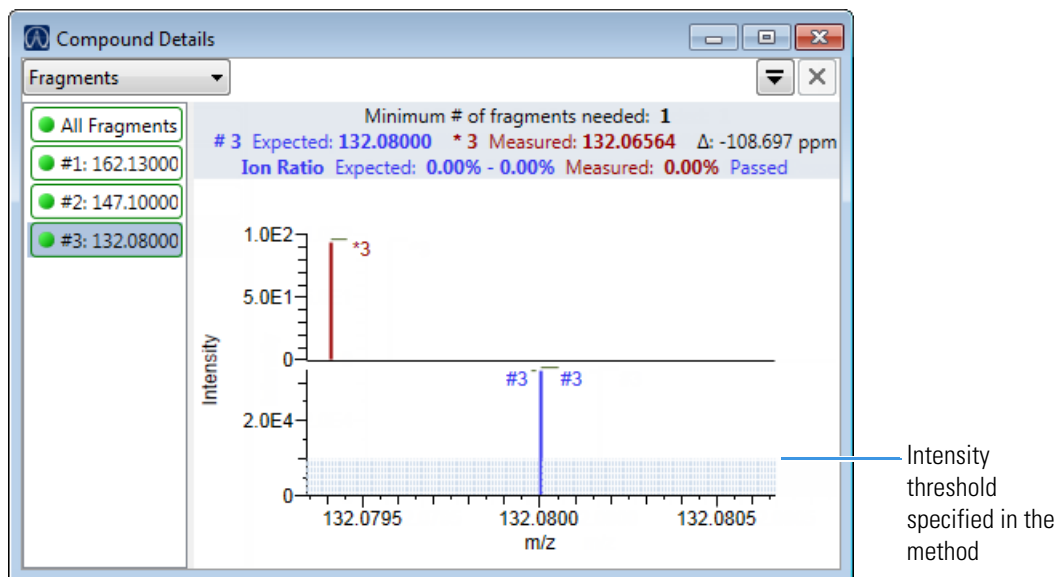


Figure 60. Fragments pane with stacked spectra for a single fragment

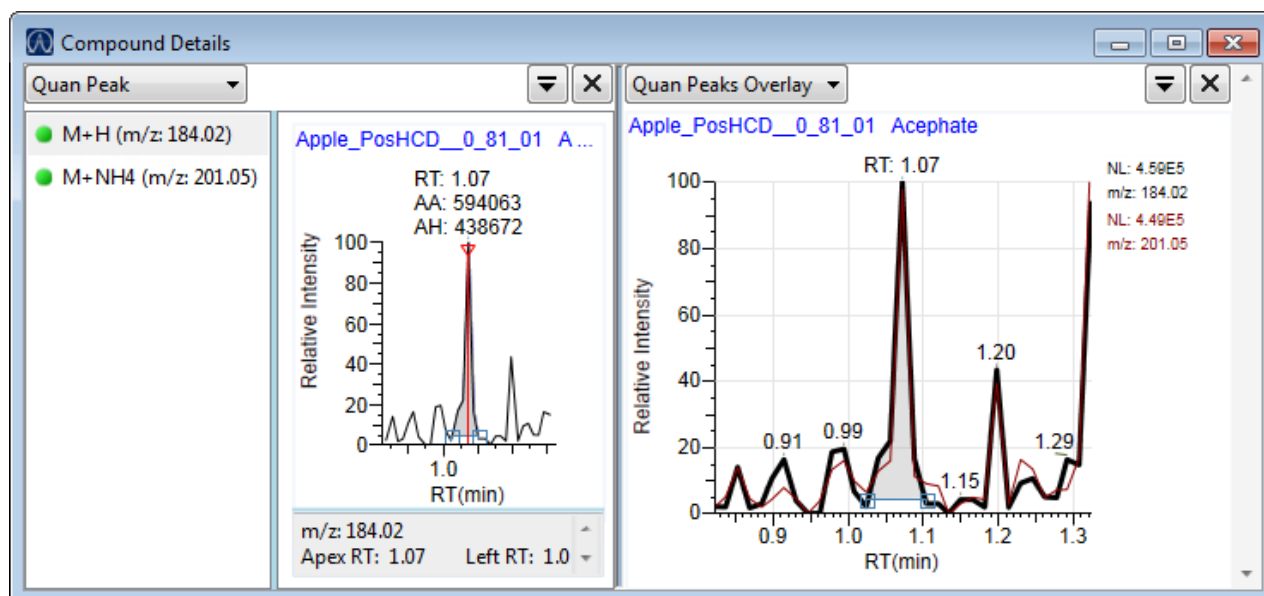


Quan Peaks Overlay

The [Quan Peaks Overlay](#) page displays overlaying quantitation peaks for compounds with multiple target peaks.

- The first target peak in the peak list appears in black in the overlay pane.
- The second target peak in the peak list appears in red in the overlay pane.
- The selected peak is bolded in the overlay pane.

Figure 61. Quan Peaks Overlay page



4 Using the Analysis Mode for Quantitation Batches

Working in Data Review for Quantitation Batches

Comparative View for Quantitation Batches

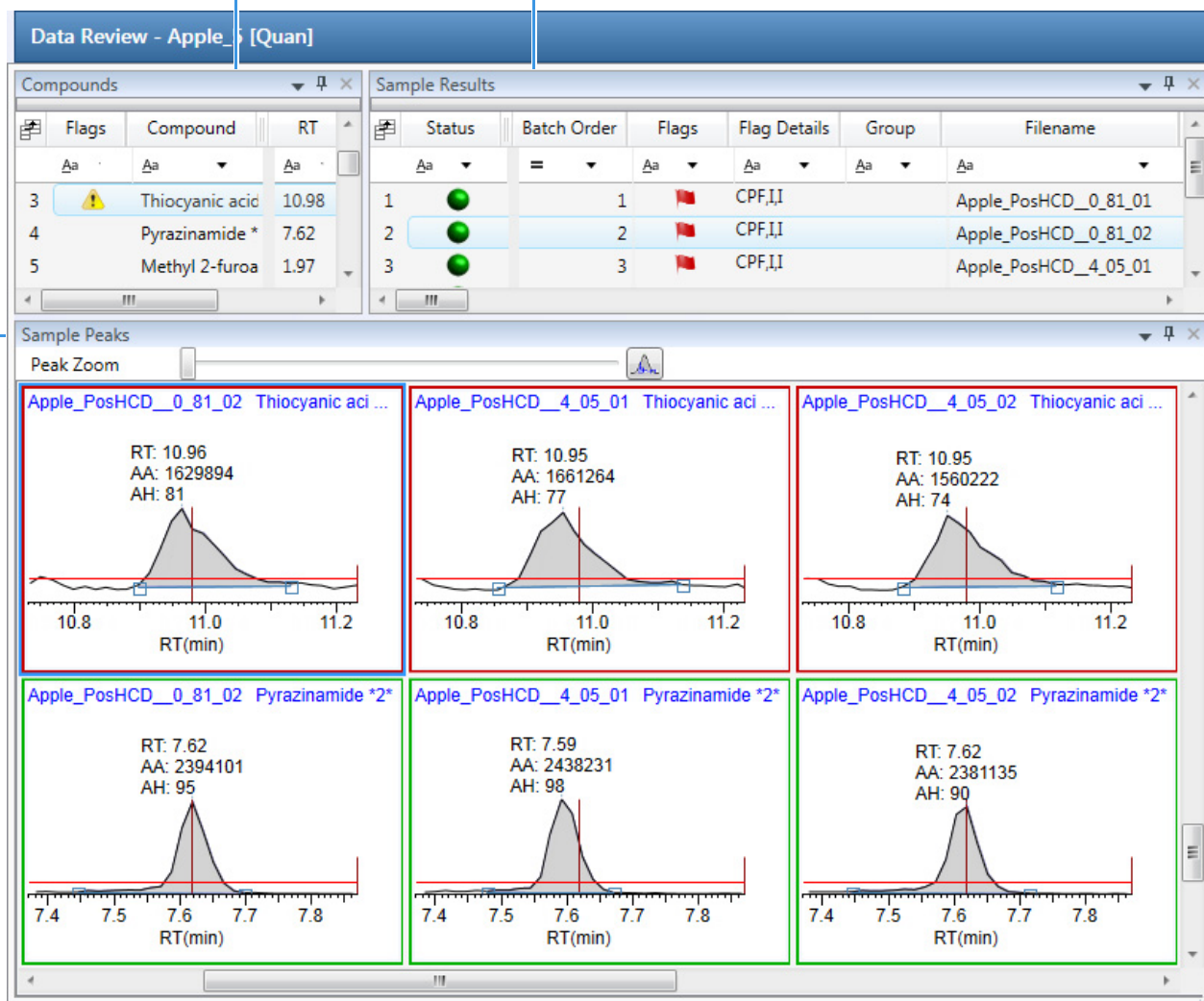
The Comparative View uses three panes to display a list of all compounds available in the method, all samples in the current batch, and the sample peak plots for all compounds found in the samples with the horizontal threshold guide.

These are the default panes and their locations:

Sample Peaks pane

Compounds pane

Sample Results pane



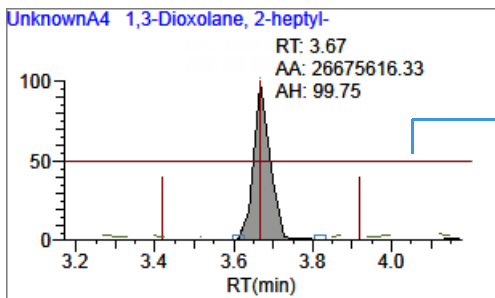
Related Topics

- [Using the Threshold Guide](#)
- [Configuring Sample Peaks Display Settings](#)
- [Manually Integrating Peaks](#)

Using the Threshold Guide

The following factors define the threshold guide that the Comparative View displays in the sample peak plots:

- The threshold method and amount that you specified in the method
- The group that you created on the Sample page
- The threshold sample that you selected on the Threshold Samples page



The horizontal guide indicates the threshold value as specified in the method.

For information about creating groups, see [Groups](#).

For information about specifying a threshold sample, see [Threshold Samples Page for Quantitation Batches](#).

For information about specifying the method to use for creating a threshold guide, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

Configuring Sample Peaks Display Settings

The Sample Peaks pane in the Comparative View displays one row per compound and one column per sample. The Sample Peaks pane displays all samples in a group when you select any of the samples belonging to that group.

For information about creating groups, see [Groups](#).

❖ To change the Sample Peaks pane display

1. From the View menu, choose **Chromatogram Pane Settings**.

The [Chromatogram Plot Settings Dialog Box](#) opens.

2. To change the number of rows or columns to fit in the Sample Peaks pane, type new values in the Number of Rows or Number of Columns boxes.

These values do not change the number of rows (compounds) and columns (samples) that are available in the Sample Peaks pane. These values determine how many rows and columns you want to view at one time in the display. The default is three rows and three columns.

4 Using the Analysis Mode for Quantitation Batches

Working in Data Review for Quantitation Batches

In the following examples, the Number of Rows and Number of Columns are set to 1 (Number of Rows equals 1 and Number of Columns equals 1) and the Number of Rows and Number of Columns are set to 4 (Number of Rows equals 4 and Number of Columns equals 4).

Figure 62. Number of Rows equals 1 and Number of Columns equals 1

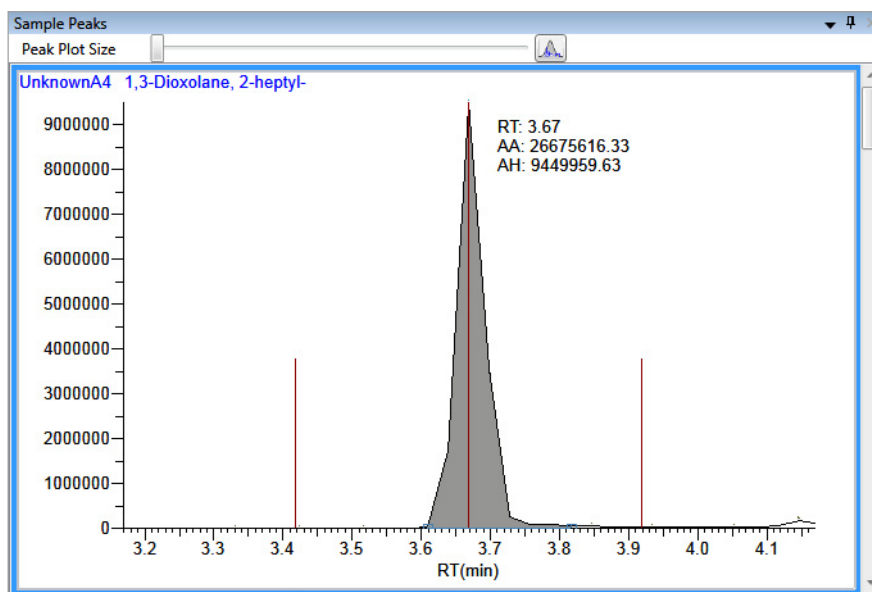
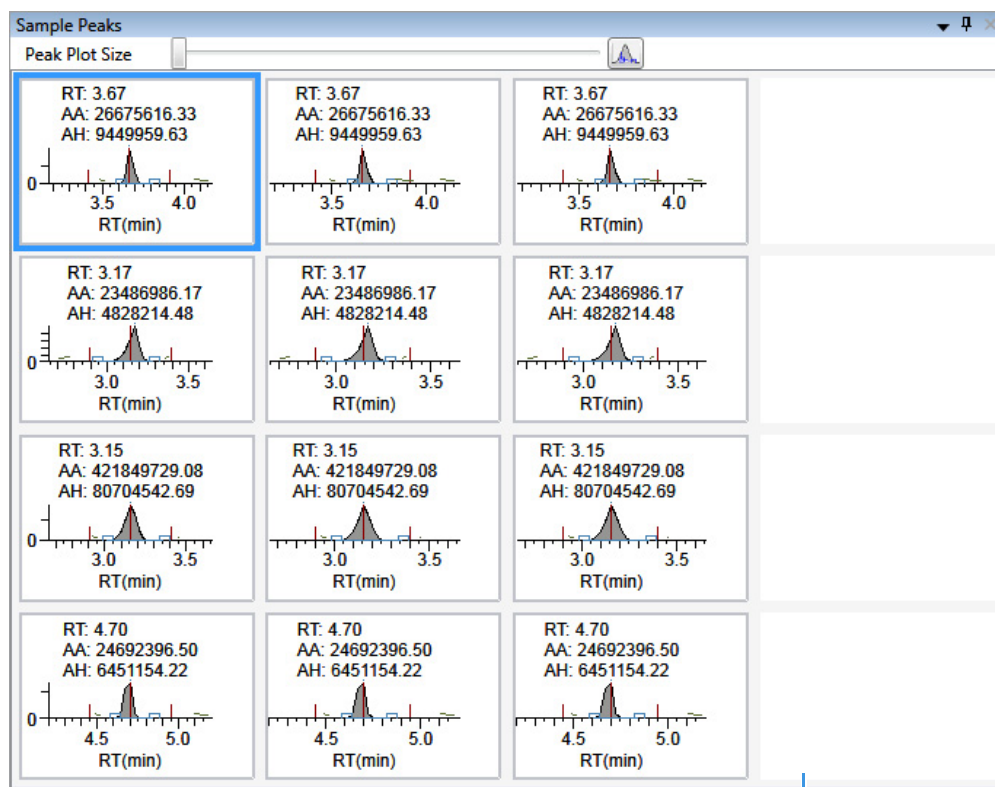


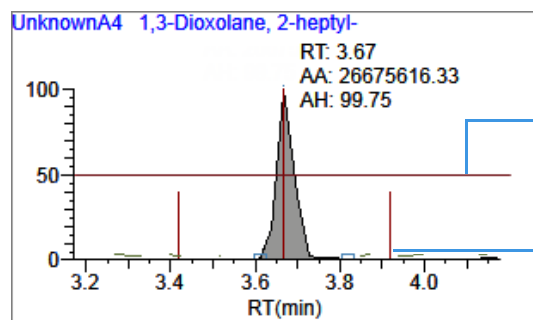
Figure 63. Number of Rows equals 4 and Number of Columns equals 4



Because there are only three samples, this column is empty.

3. To change the display type for the y -axis scale, select one of the following:

- **Relative:** Displays the y -axis scale from 0 through 100.



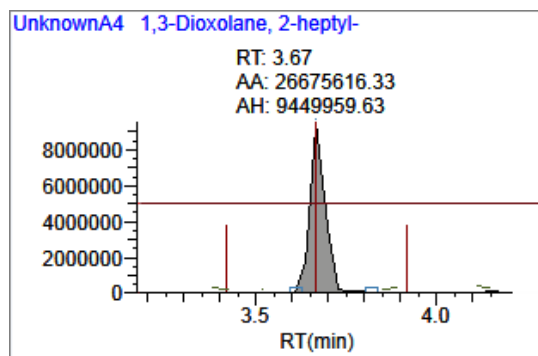
The horizontal bar indicates the threshold value as specified in the method.

Vertical bars indicate the expected retention time and window, as specified in the method.

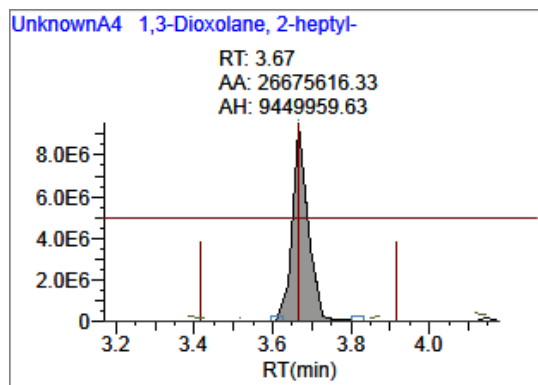
4 Using the Analysis Mode for Quantitation Batches

Working in Data Review for Quantitation Batches

- **Absolute:** Displays the y -axis scale from 0 to the actual value of the most intense peak in the group.



- **Label in Scientific Notation:** Displays the y -axis scale in scientific notation.



Note The Sample Peaks pane displays a y axis only in the first chromatogram in each row. The limits of the scale are determined by the most intense peak in the group.

4. Specify which labels you want to display in the sample peak plots.

For an example of all available peak plot labels, see [Peak Plot Labels](#).

Chromatogram Plot Settings Dialog Box

Use the Chromatogram Plot Settings dialog box to change the Sample Peaks pane display.

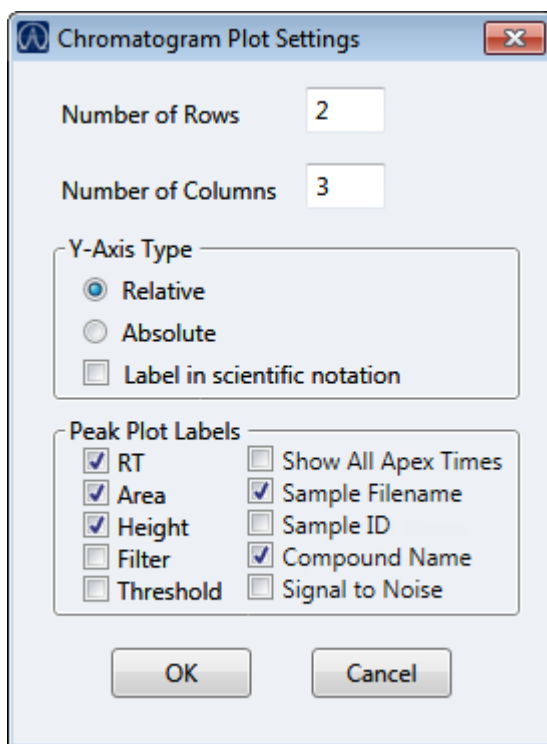


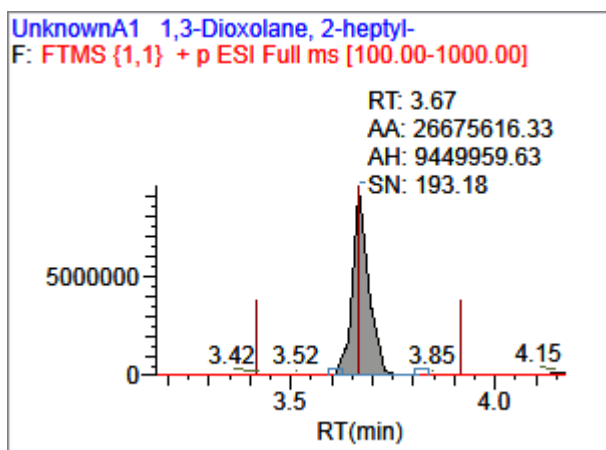
Table 35. Chromatogram Plot Settings dialog box parameters (Sheet 1 of 2)

Parameter	Description
Number of Rows	Specifies the number of rows visible in the Sample Peaks pane. When Number of Rows equals 1, the application scales the height of all chromatograms to fill the Y dimension of the Sample Peaks pane. Default: 3
Number of Columns	Specifies the number of columns visible in the Sample Peaks pane. When Number of Columns equals 1, the application scales the width of all chromatograms to fill the X dimension of the Sample Peaks pane. Default: 3
Y-Axis Type	Displays the y-axis scale as Relative (to the most intense peak), Absolute, or in scientific notation.

Table 35. Chromatogram Plot Settings dialog box parameters (Sheet 2 of 2)

Parameter	Description
Peak Plot Labels	Displays or hides the following peak labels: <ul style="list-style-type: none"> • RT • Area • Height • Filter • Threshold • Show All Apex Times • Sample Filename • Sample ID • Compound Name • Signal to Noise

Example with all peak labels displayed:

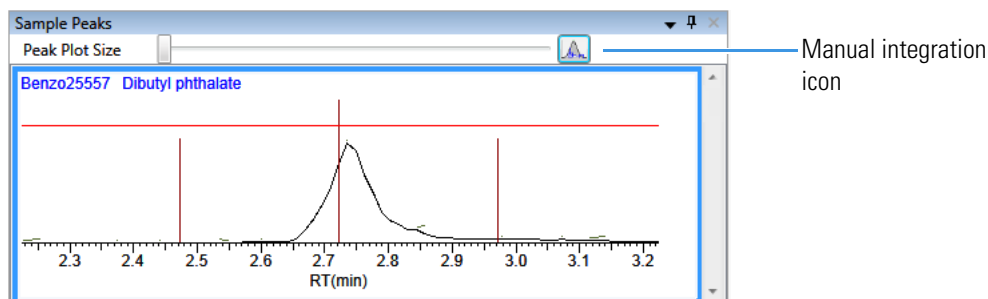


Manually Integrating Peaks

Use the manual integration feature to manually add a peak. You can manually add a peak in a chromatogram plot only when the application fails to identify a peak.

❖ To manually integrate a peak

1. In the Sample Peaks plot, click the **Manual Integration** icon.



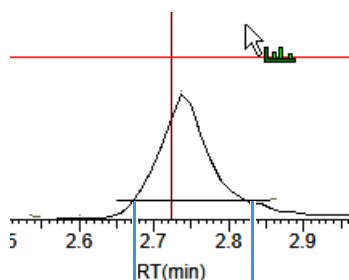
The cursor changes to look like this:



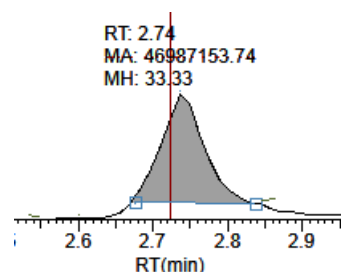
2. Integrate a peak as follows:
 - a. Drag the cursor to define the beginning and ending base points for the new peak.

Note You must drag the cursor inside the *x* axis and *y* axis.

- b. Click outside the plot to refresh the display.



Click the first base point, and then drag to the second base point.



Click outside the plot to refresh the view.

The application identifies the peak and indicates the manual integration in the labels.

3. To zoom in on an area, do the following:

- a. Drag the cursor below the *x* axis or to the left of the *y* axis.

The plot zooms to fit the defined X or Y dimension into the entire pane. The application zooms all compounds in the row to the same scale.

- b. To return to the original view, right-click and choose **Reset Scaling**.

Qualitative View for Quantitation Batches

The Qualitative View uses several different panes to display qualitative information for the selected sample. See [Displaying Qualitative View Panes](#).

If the application finds no detected peaks for the selected sample, you can manually add peaks.

To see processed data for a sample in the Qualitative View, you must select the Qual Processing parameter for that sample in the Batch View before you process the batch. See [Batch View Sample List](#).

These are the default panes and their locations:

The Qualitative View displays data in the following panes:

- [Samples Pane](#)
- [Peaks Pane](#)
- [Sample Chromatogram Pane](#)
- [Peak Chromatogram Pane](#)
- [Spectrum Pane \(Library and Data\)](#)
- [Library Hits Pane](#)

Displaying Qualitative View Panes

The Qualitative View displays data using multiple panes: Samples, Peaks, Sample Chromatogram, Peak Chromatogram, Spectrum, and Library Hits. You can display, hide, or move any of these panes.

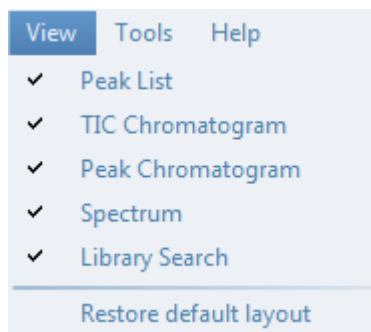
❖ To display or hide a Qualitative View pane

From the View menu, choose to display or hide the following:

- **Peak List:** Displays the Peaks pane
- **TIC Chromatogram:** Displays the Sample Chromatogram pane
- **Peak Chromatogram:** Displays the Peak Chromatogram pane
- **Spectrum:** Displays the Spectrum pane
- **Library Search:** Displays the Library Hits pane

Note The Samples pane is required for the Qualitative View display. You cannot hide the Samples pane.

A check mark indicates a displayed pane.




For procedures about creating docked, floating, or tabbed panes, see [Moving Data Review Panes](#).

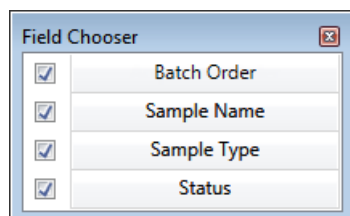
Samples Pane

Use the Samples pane in the Qualitative View to select a specific sample. The associated [Peaks Pane](#) displays all peaks found in the selected sample.

❖ To hide or display columns in the Samples pane

1. Click the **Field Chooser** icon, , in the upper left corner of the pane.

The Field Chooser displays all available columns of data for the Samples pane.



2. Select the check box for each column that you want to display, or clear the check box for each column that you want to hide.

The application immediately displays or hides the column in the Samples pane.


3. Click  to close the Field Chooser.

Figure 64. Samples pane







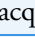

Samples				
	Batch Order	Status	Sample Name	Sample Type
= ▼	Aa	Aa	Aa ▼	Aa ▼
1	1	●	Apple_PosHCD_0_81_01	Specimen
2	2	●	Apple_PosHCD_0_81_02	Specimen
3	3	●	Apple_PosHCD_4_05_01	Specimen

Table 36. Samples pane columns

Column	Description
Batch Order	Sequentially numbers the samples.
Status	 Sample is not acquired.  Sample is acquired but not processed.  Sample is acquired and processed.  Sample is currently acquiring. <p>Note When you include unknown screening features in the quantitation method and you choose to process with only the quantitation criteria, the Sample View shows the Status for the samples as acquired and processed () , and the Unknown Screening View shows the Status for the samples as acquired but not processed () .</p> <p>Note When you include unknown screening features in the quantitation method and you choose to process with only the unknown screening criteria, both the Sample View and the Unknown Screening View show the Status for the samples as acquired and processed () .</p>
Sample Name	A user-defined name that identifies a sample.
Sample Type	Defines how the application processes the sample data. Each sample is classified as one of the following sample types: Solvent, Specimen, QC, Calibrator, Hydrolysis, Unextracted, or Negative.

Peaks Pane

The [Peaks Pane](#) in the Qualitative View works with the Samples pane to display graphical values for a unique sample and peak combination.

Follow these procedures:

- [To display peaks for a specific compound](#)
- [To hide or display columns in the Peaks pane](#)

❖ To display peaks for a specific compound

1. From the Samples pane, select a sample.

The Peaks pane displays the retention times for peaks identified in the selected sample, the values for the best match methods for each peak, and the library match.

The method specifies which technique to use for identifying peaks: peaks within a specific retention time range, as a minimum percentage of the height or area of the largest peak, or as a minimum percentage of the nearest internal standard peak. You can change the method for identifying peaks in the Method Template Editor. Refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

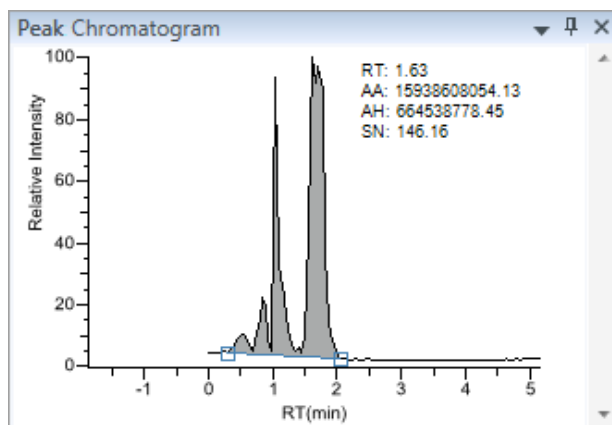
2. In the Peaks pane, select a peak in the sample.

Peaks						
Filter: + c ESI Full ms [200.00-800.00]						
Peak RT	SI	RSI	MP	Est Amt	Library Hit	
9.82	124	732	0	0.000	Phenomorphane	
1.63	161	570	4	0.000	Sulfasomizole	

The application displays the selected peak in the [Peak Chromatogram](#) pane, displays the Qual Data and Qual Library sections in the [Spectrum](#) pane, and locates the selected peak in the [Sample Chromatogram](#) pane.

- The Qual Data section shows spectrum data for the peak in the raw data file.
- The Qual Library section shows actual spectrum for the identified library compound.

Figure 65. Peak Chromatogram pane



Note When you select a data-dependent sample, the peak can be from either a full scan or a QED spectrum of an SRM-filtered chromatogram.

Figure 66. Spectrum pane

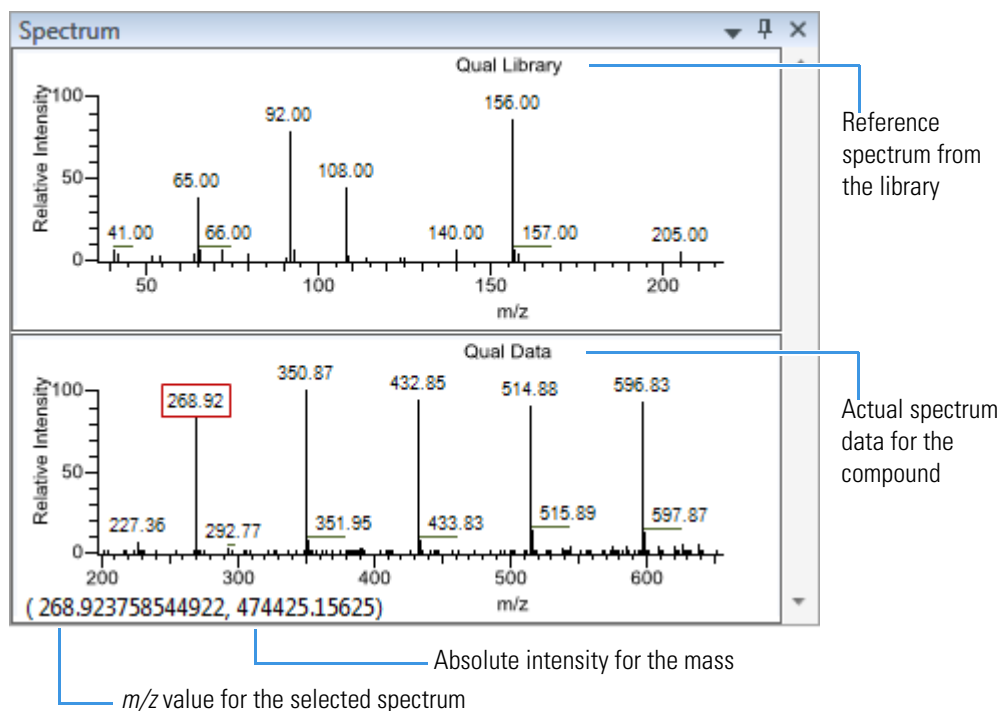
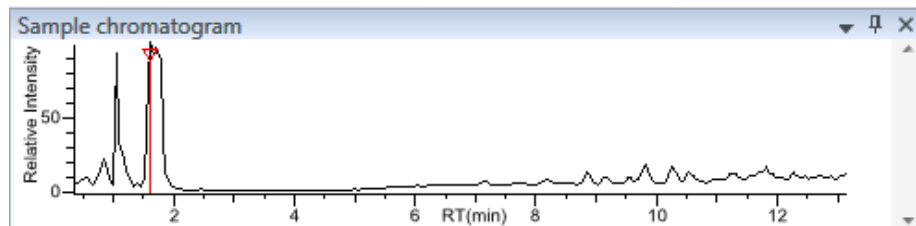



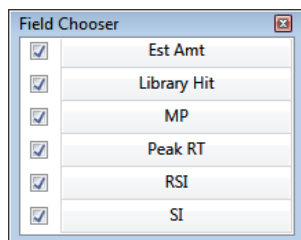
Figure 67. Selected peak in the Sample Chromatogram pane



❖ To hide or display columns in the Peaks pane


1. Click the **Field Chooser** icon, , in the upper left corner of the pane.

The Field Chooser displays all available columns of data for the Peaks pane.



2. Select the check box for each column that you want to display, or clear the check box for each column that you want to hide.

The application immediately displays or hides the column in the Peaks pane.

3. When you are finished modifying the column display, click  to close the Field Chooser.

Peaks Pane

Use the features in the Peaks pane to display graphical values for each sample and peak combination.

Figure 68. Peaks pane

Peak RT	SI	RSI	MP	Est Amt	Library Hit
9.82	124	732	0	0.000	Phenomorphane
1.63	161	570	4	0.000	Sulfasomizole

Table 37. Peaks pane parameters

Parameter	Description
Filter	Filter used to identify the peaks. Specified in the raw data file or the master method. When your raw data file is data-dependent, the filter indicates this with a “d”. <div> <div>Filter: + c d ESI Full ms [200.00-800.00]</div> <div>Data-dependent filter</div> </div>
Peak RT	Peak retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column.
SI	Search index method used to search the NIST library.
RSI	(Reverse search index) A method used to search the NIST library. A reverse search compares a library entry to an unknown compound (whereas a forward search compares the mass spectrum of an unknown compound to a mass spectral library entry).
MP	Match probability.
Est Amt	Estimated amount of the compound.
Library Hit	Library compound that matches the identified peak.

Sample Chromatogram Pane

The [Sample Chromatogram Pane](#) in the Qualitative View displays all peaks in the selected sample. The peak selected in the Peaks pane displays a red marker.

❖ To zoom in on a peak

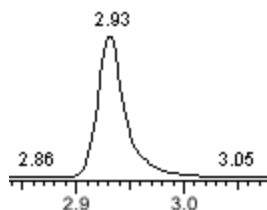
1. In the Sample Chromatogram pane, drag the cursor to delineate a rectangle around the peak.

The delineated area expands to fill the view.

2. To restore the default view, right-click the Sample Chromatogram pane and choose **Reset Scaling**.

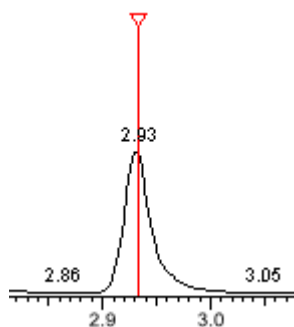
❖ To manually add a peak

1. Zoom in to better identify which peak to add to the results set.



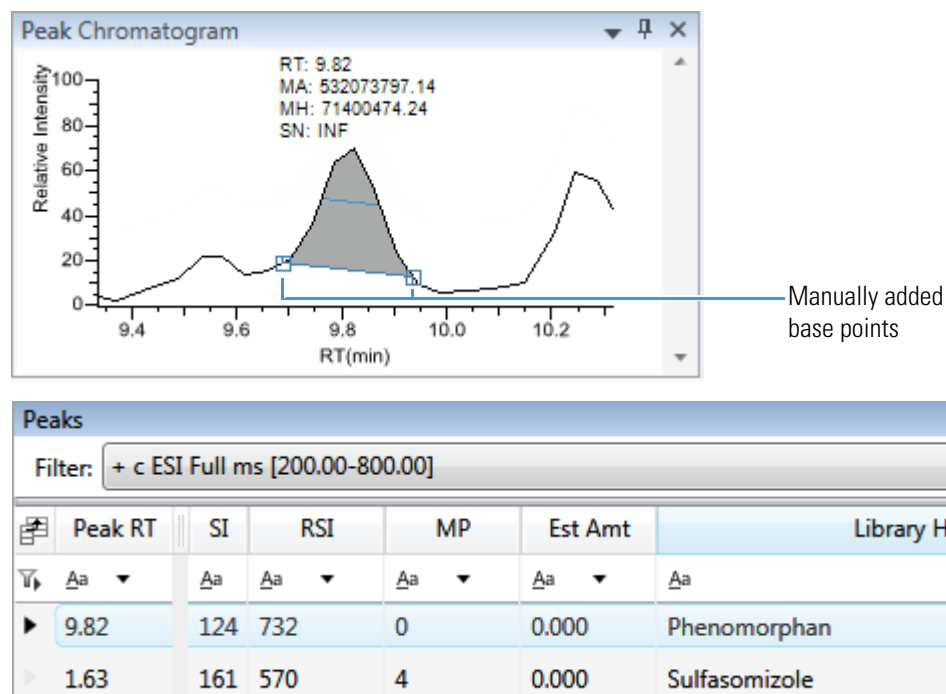
2. Right-click the Sample Chromatogram pane, and choose **Add Qual Peak**.
3. Click to indicate the left and right base points for the peak.

The application marks the peak in the Sample Chromatogram pane.



The application places the peak delimiter tags at the base point locations and automatically updates the peak values in the Peaks pane and Peak Chromatogram pane. See [Peak Chromatogram pane with a manually added peak](#).

Figure 69. Peak Chromatogram pane with a manually added peak



Sample Chromatogram Pane

Use the features in the Sample Chromatogram pane to display peaks in the selected sample.

Figure 70. Sample Chromatogram pane

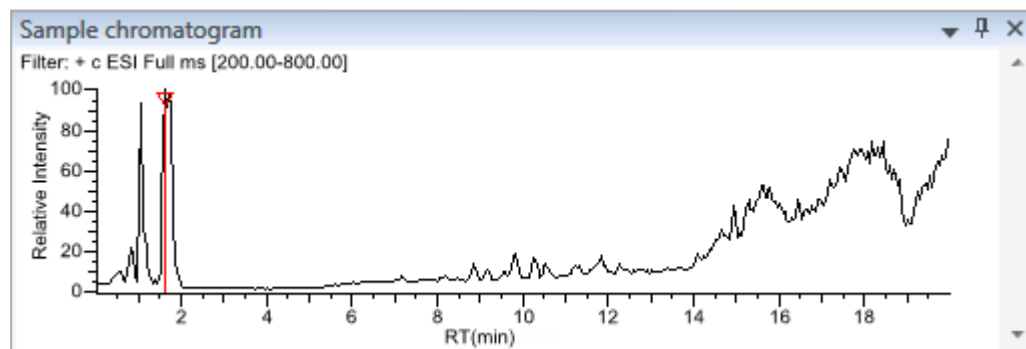


Table 38. Sample chromatogram pane shortcut menu commands

Command	Description
Add Qual Peak	Select the beginning and ending base points for a new qualitative peak. Available only when no peak is detected.
Reset Scaling	Resets the original scaling after a zoom operation.
Copy to Clipboard	Copies the graphic display to the Clipboard.

Peak Chromatogram Pane

The [Peak Chromatogram Pane](#) in the Qualitative View displays the selected peak.

Follow these procedures:

- [To zoom in on a peak](#)
- [To remove a peak](#)
- [To switch between method and manual integration modes](#)
- [To customize the labels in the display](#)

❖ To zoom in on a peak

1. In the chromatogram plot, drag the cursor to delineate a rectangle around the peak.

The delineated area expands to fill the view.

2. To restore the default view, right-click the chromatogram plot and choose **Reset Scaling**.

❖ To remove a peak

Right-click the Peak Chromatogram pane, and choose **Remove Qual Peak**.

The application removes the peak displayed in the Peak Chromatogram pane. All data for this peak are removed from the Qualitative View panes.

❖ To switch between method and manual integration modes

Right-click the Peak Chromatogram pane and choose either **Method Integration** or **Manual Integration**.

Initially, the method integration and manual integration settings that are stored for a compound and file are identical. When you switch modes, the saved result set does not change. However, when manual data are available, both the Peak Chromatogram plots and the result set update as you switch between method and manual modes.

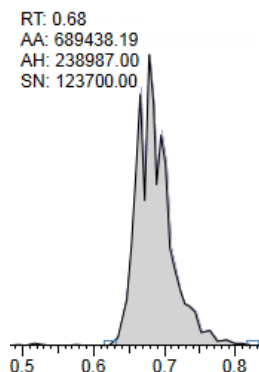
As you switch between modes, each pane reflects the changes. The generated reports for the data identify the manual modifications.

❖ To customize the labels in the display

1. Right-click and hold the cursor over **Peak Labels** in the shortcut menu.
2. To display labels for the peak retention time (RT), area (AA), height (AH), or signal-to-noise (SN), select the appropriate check box.

See [Example of displayed labels](#).

Figure 71. Example of displayed labels



Peak Chromatogram Pane

Use the features in the Peak Chromatogram pane to display the selected peak.

Figure 72. Peak Chromatogram pane

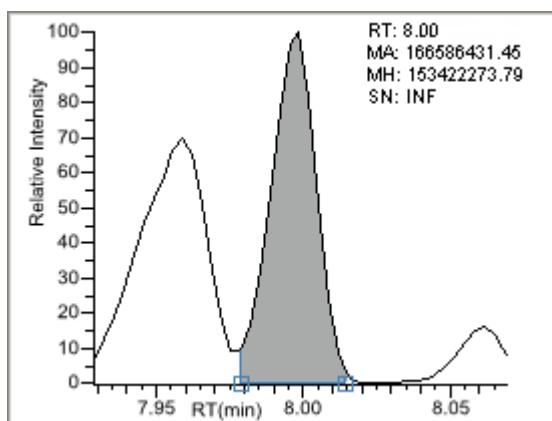


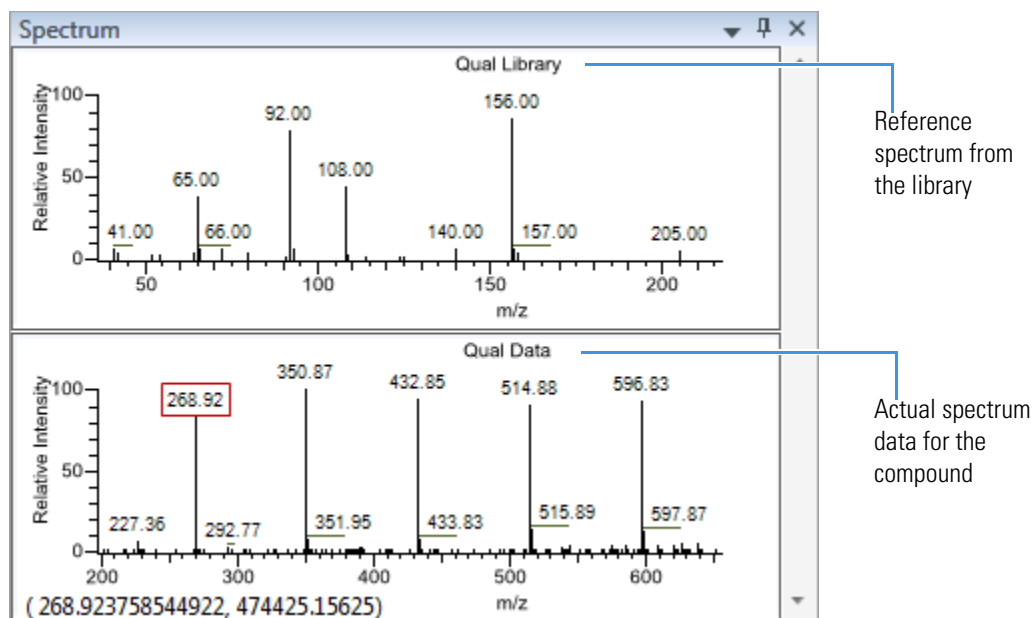
Table 39. Peak Chromatogram pane shortcut menu commands

Command	Description
Reset Scaling	Resets the original scaling after a zoom operation.
Method Integration	Displays method integration settings.
Manual Integration	Displays manual integration settings.
Peak Labels	Displays or hides the peak labels (Label Area, Label Retention Time, Label Height, or Label to Noise).
Remove Qual Peak	Available only for manually added peaks. Removes the peak displayed in the Peak Chromatogram pane.
Copy to Clipboard	Copies the graphic display to the Clipboard.

Spectrum Pane (Library and Data)

The Spectrum pane in the Qualitative View displays the reference spectrum from the library and the spectrum data for the selected sample. The top pane displays the spectrum for the identified compound found in the reference library; the bottom pane displays the actual spectrum data for the selected peak.

Figure 73. Spectrum pane



❖ To zoom in on a peak

1. In either spectrum plot, drag the cursor to delineate a rectangle around the peak.

The delineated area expands to fill the view.

2. To restore the default view, right-click the spectrum plot and choose **Reset Scaling**.

❖ To display detailed m/z values

In the spectrum pane, hold the cursor over the m/z value for a peak.

The application indicates the rounded-off m/z value with a red box and displays the complete value in the lower left corner of the display.

Library Hits Pane

The **Library Hits pane** in the Qualitative View displays the best library matches for the selected peak. Use this pane to select a different library entry for the peak.

When you select a library entry other than the original entry, the TIC Report and TIC Summary Report indicate this with a “P” flag:

Peak:	Retention Time	Area	Height	Inj Estimate	In-sample Est	Flag
Naphthalene	7.95	13829174	10605061	0.000	0.000	P

P flag

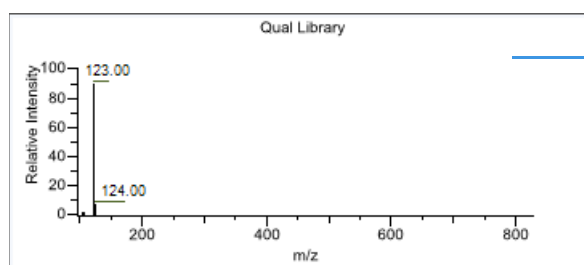
❖ To change the library entry for a selected peak

In the Library Hits pane, select the check box for the library entry that you want to use to identify the selected peak.

- In the Spectrum pane, the reference spectra change to show the spectra for the selected library entry.
- In the Peaks pane, the SI, RSI, MP, and Compound values update to reflect the selected library entry.

Library hits					
	Rank	SI	RSI	MP	Library entry
<input type="radio"/>		=	=	=	Aa
<input checked="" type="radio"/>	1	332	978	0	2-Hexanone
<input type="radio"/>	2	320	966	0	Succinic anhydride
<input type="radio"/>	3	314	959	0	Propane, 1-(ethenylloxy)-

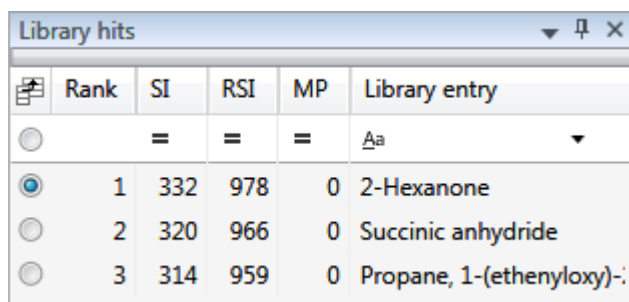
Selected library entry
in the Library Hits pane



Reference spectra
for Hexanone

	Peak RT	SI	RSI	MP	Est Amt	Library Hit
	Aa	Aa	Aa	Aa	Aa	Aa
▶	0.00	332	978	0	0.000	2-Hexanone
▶	4.46	145	860	0	0.000	Methane, bis(2-chloroethoxy)-

Peak list for
Hexanone

Figure 74. Library Hits pane


<input type="checkbox"/>	Rank	SI	RSI	MP	Library entry
<input type="radio"/>	=	=	=	=	Aa
<input checked="" type="radio"/>	1	332	978	0	2-Hexanone
<input type="radio"/>	2	320	966	0	Succinic anhydride
<input type="radio"/>	3	314	959	0	Propane, 1-(ethenyloxy)-

Table 40. Library Hits pane parameters

Parameter	Description
<Check box column>	Indicates selected library entries for the selected peak.
Rank	Indicates the order of best matches between the selected peak and library entries.
SI	(Search index) A method used to search the NIST library.
RSI	(Reverse search index) A method used to search the NIST library. A reverse search compares a library entry to an unknown compound (whereas a forward search compares the mass spectrum of an unknown compound to a mass spectral library entry).
MP	Match probability.
Library Entry	Library compound that matches the identified peak.

Column Parameters for Compound Results and Sample Results

The Compound Results and Sample Results data review pages use the following parameters and features:

- [Common Column Parameters](#)
- [Inactive and Excluded Compounds](#)
- [Exporting Compounds](#)

Common Column Parameters

Table 41. Common parameters for Compound Results and Sample Results tables (Sheet 1 of 8)

Column	Description
%CV	Coefficient of Variance. Standard deviation of the multiple samples of one level, multiplied by 100, and divided by the average of the multiple samples of that level. This calculation is based on the areas of the peaks.
%Diff	The calculated amount minus the expected amount, divided by the expected amount, and then multiplied by 100.
%RSD	Standard deviation of the multiple samples of one level, multiplied by 100, and divided by the average of the multiple samples of that level. This calculation is based on the calculated amounts. Note This RSD value is not the same as the RSD value used with the Average RF curve type in the method. Refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the <i>TraceFinder Lab Director User Guide</i> . The application uses this %RSD value in Data Review and in the Compound Calibration Report when you acquire multiple samples for the same QC or Calibrator samples.
Active	Displays or hides a compound for a particular sample. <ul style="list-style-type: none"> • When a compound is marked inactive, the application does not remove its data and calculated values from the result set. Instead, the application masks the appearance of that compound for that particular sample and grays the compound in the compounds list. • When a calibration standard is marked inactive, the application no longer uses the data file’s calibration point for the calibration and removes it from the graphical view of the calibration curve displayed in the Compound Details pane. It is no longer part of the result set. <p>In a Sample View, the Active parameter is in the Compound Results pane.</p> <p>In a Compound View, the Active parameter is in the Sample Results pane.</p>
Actual RT	Actual retention time for the compound. Retention time is the time after injection when a compound elutes and the total time that the compound is retained on the chromatograph column.

Table 41. Common parameters for Compound Results and Sample Results tables (Sheet 2 of 8)

Column	Description
Adduct	The most intense adduct for the retention time for a compound.
Area	The area obtained by integrating peak intensities from the start to the end of the peak. When the Response Ratio is specified as Area, this column displays an asterisk (*Area).
Batch Order (Sample Results only)	Sequentially numbers the samples in the batch.
Calculated Amt	The amount present in the sample, as determined using the calibration curve and the response ratio.
Channel	Specifies the channel on which the sample was run. If the sample is not acquired, the value is Pending. The Channel column is available only when you have activated multiplexing in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the <i>TraceFinder Lab Director User Guide</i> .
Compound (Compound Results only)	Name of the identified compound.
Comments (Sample Results only)	A user-defined comment for the sample.
Confirm	Specifies that the compound meets all required processing criteria. Refer to the instructions “To display only confirmed peaks” in Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the <i>TraceFinder Lab Director User Guide</i> .

To display the confirmation criteria, hold your cursor over the indicator.

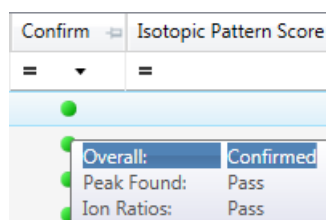


Table 41. Common parameters for Compound Results and Sample Results tables (Sheet 3 of 8)

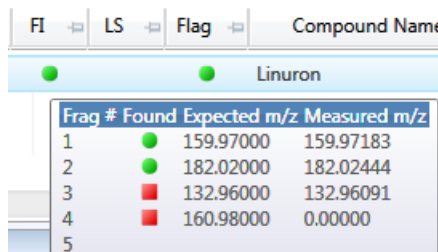
Column	Description																								
FI	<p>Fragment Ions flag. The application displays one of these indicators:</p> <ul style="list-style-type: none">A green circle (pass) when the measured <i>m/z</i> value of any of the fragments is within the mass tolerance specified in the method. On the Isotopes page in the Spectrum pane, the All Isotopes and Multi-Isotopes flags are also green.A red square (fail) when the measured <i>m/z</i> value of none of the fragments is within the mass tolerance specified in the method. On the Isotopes page in the Spectrum pane, the All Isotopes and Multi-Isotopes flags are also red.A blank when there are no fragments detected. <p>To display a list of fragments and their pass/fail status, hold your cursor over the indicator.</p>  <p>The screenshot shows a table for the compound Linuron. The table has columns: Frag #, Found, Expected <i>m/z</i>, and Measured <i>m/z</i>. The data is as follows:</p> <table><thead><tr><th>Frag #</th><th>Found</th><th>Expected <i>m/z</i></th><th>Measured <i>m/z</i></th></tr></thead><tbody><tr><td>1</td><td>Green circle</td><td>159.97000</td><td>159.97183</td></tr><tr><td>2</td><td>Green circle</td><td>182.02000</td><td>182.02444</td></tr><tr><td>3</td><td>Red square</td><td>132.96000</td><td>132.96091</td></tr><tr><td>4</td><td>Red square</td><td>160.98000</td><td>0.00000</td></tr><tr><td>5</td><td></td><td></td><td></td></tr></tbody></table>	Frag #	Found	Expected <i>m/z</i>	Measured <i>m/z</i>	1	Green circle	159.97000	159.97183	2	Green circle	182.02000	182.02444	3	Red square	132.96000	132.96091	4	Red square	160.98000	0.00000	5			
Frag #	Found	Expected <i>m/z</i>	Measured <i>m/z</i>																						
1	Green circle	159.97000	159.97183																						
2	Green circle	182.02000	182.02444																						
3	Red square	132.96000	132.96091																						
4	Red square	160.98000	0.00000																						
5																									
Filename	A user-defined name that identifies a sample.																								
Final Units	<p>Specifies the calculated amount.</p> <p>Default: 1</p>																								
Flag Details	<p>Indicates all errors found in the compound.</p> <p>Type of error:</p> <ul style="list-style-type: none">I: Confirming ion coelution failureA: Amount errorB: Matrix blank errorPK: Peak not foundLS: Library matching errorIP: Isotope errorFI: Fragment ions errorIR: Ion ratio error																								
Flags	Caution flag displayed when a compound within the sample has an error.																								
Formula	The formula for the peak as specified in the compound database.																								
Group (Sample Results only)	Threshold group to which a sample belongs. You can view samples by group in the Comparative View.																								
Group Average (Compound Results only)	Average of the samples in a threshold group. You can view averaged samples in the Compound View.																								

Table 41. Common parameters for Compound Results and Sample Results tables (Sheet 4 of 8)

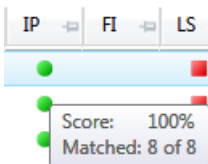
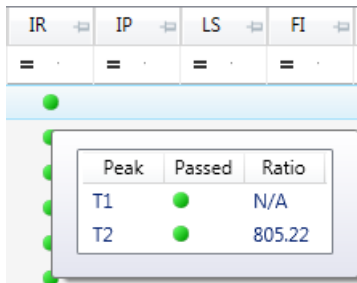
Column	Description
Height	Specifies the distance from the peak maximum to the peak base, measured perpendicular to the ordinate. When the Response Ratio is specified as Height, this column displays an asterisk (*Height).
Integration Mode	Specifies the integration mode (Method, Manual, or User). See Quan Peak .
Ion Ratio	Specifies the ratio of the confirming ion's response to the target ion's response.
Ion Type	Specified as b or y with a numerical designation. Available only for peptide databases.
IP	<p>Isotopic Pattern flag. The application displays one of these indicators:</p> <ul style="list-style-type: none"> • A green circle (pass) when the score percentage is higher than the specified fit threshold percentage • A red square (fail) when the score percentage is lower than the specified fit threshold percentage • A blank when the parameter is not scored <p>To display the score of matched isotopes, hold your cursor over the indicator.</p> 
IR	<p>Ion Ratio flag. The application displays one of these indicators:</p> <ul style="list-style-type: none"> • A green circle when the ion ratio is within the acceptable ion ratio range • A red square when the ion ratio is not within the acceptable ion ratio range <p>To display the ion ratio results, hold your cursor over the indicator.</p> 
Isotopic Pattern Score (%)	The percentage of the number of total isotopes to the number of matched isotopes.
ISTD Amt	Amount of internal standard.
ISTD Response	Response of the internal standard.
Level (Sample Results only)	The level defined for a calibration sample or quality control sample.

Table 41. Common parameters for Compound Results and Sample Results tables (Sheet 5 of 8)

Column	Description
Lib Match Name	<p>The name of the best matching compound in the library search. When the application finds a match in the library, this column displays the matching library entry with the highest score.</p> <ul style="list-style-type: none"> When the application does not perform a library search, this column displays “N/A” in black text. When the application does not perform an MS/MS scan, this column displays “N/A” in red text.
Library Match Rank	<p>Displays the ranking of the library match. When the application finds a match in the library, this column displays the library entry’s relative rank, in the format “x of y”, where</p> <ul style="list-style-type: none"> x = the rank of the highest scoring library match. y = the total number of library matches from the list of matches for a particular adduct that contains the highest scoring match. <p>Results are as follows when the application performs both a library search and an MS/MS scan and both the library entry and the formula match the target compound:</p> <ul style="list-style-type: none"> A library score that is higher than or equal to the score threshold meets the criteria, and the values in this column appear in green. A library score that is lower than the score threshold fails to meet the criteria, and the values in this column appear in red. <p>When the application does not perform a library search, this column displays “N/A” in black text.</p> <p>When the application does not perform an MS/MS scan, this column displays “N/A” in red text.</p>
Library Probability Percentage	Specifies the probability of the match in the NIST library for forward searches only.

Table 41. Common parameters for Compound Results and Sample Results tables (Sheet 6 of 8)

Column	Description
Library Score (%)	<p>Specifies the score from the library fit. When the application finds a match in the high resolution NIST library or in the mzVault library, this column displays the highest score associated with the Lib Match Name parameter.</p> <p>Results are as follows when the application performs both a library search and an MS/MS scan and both the library entry and the formula match the target compound:</p> <ul style="list-style-type: none"> • A library score that is higher than or equal to the score threshold meets the criteria, and the values in this column appear in green. • A library score that is lower than the score threshold fails to meet the criteria, and the values in this column appear in red. <p>When the application does not perform a library search, this column displays “N/A” in black text.</p> <p>When the application does not perform an MS/MS scan, this column displays “N/A” in red text.</p> <p>Valid range: 1 through 100%</p>
LS	<p>Library Search flag. The application displays one of these flags:</p> <ul style="list-style-type: none"> • A green circle when the library search is successful. • A red square when the library search is not successful.
<i>m/z</i> (Apex)	<p>Mass-to-charge ratio found in the spectra for the peak. Assumes that the charge is 1.</p> <p>When the application successfully integrates the peak, this column displays the charged <i>m/z</i> value for the compound, which is the highest intensity in the apex scan.</p> <p>When the application cannot successfully integrate the peak, this column displays “N/F” (not found).</p>
<i>m/z</i> (Delta)	<p>Difference between the <i>m/z</i> (Expected) and <i>m/z</i> (Apex). Assumes that the charge is 1.</p> <p>When the <i>m/z</i> (Apex) column displays <i>m/z</i> value for the compound, this column displays the delta <i>m/z</i> corresponding to the highest intensity in the apex scan.</p> <ul style="list-style-type: none"> • When the mass tolerance is specified in ppm in the master method, then $m/z \text{ (Delta)} = 1\,000\,000 \times ([m/z \text{ (Apex)} - m/z \text{ (Expected)}] \div m/z \text{ (Expected)})$. • When the mass tolerance is specified in mmu in the master method, then $m/z \text{ (Delta)} = 1000 \times (m/z \text{ (Apex)} - m/z \text{ (Expected)})$.

Table 41. Common parameters for Compound Results and Sample Results tables (Sheet 7 of 8)

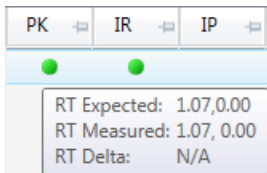
Column	Description
<i>m/z</i> (Expected)	<p>Mass-to-charge ratio from the compound database. Assumes that the charge is 1.</p> <ul style="list-style-type: none"> When an adduct is found, the application displays the neutral mass value for the compound (calculated from the neutral formula) ± the mass of the most intense adduct ion found for the compound. When no adduct is found, the application displays the neutral mass value for the compound ± the mass of the first adduct entered in the compound database. <p>For details about defining adducts for the compound database, refer to Chapter 1, “Using the Configuration Console,” in the <i>TraceFinder Lab Director User Guide</i>.</p> <p>For details about adding adducts to compounds, refer to Chapter 2, “Using Compound Databases in the Method Development Mode,” in the <i>TraceFinder Lab Director User Guide</i>.</p> <p>Note When the adduct is a gain, the adduct mass is a positive number. When the adduct is a loss, the adduct mass is a negative number. The resulting mass value after adding or subtracting the adduct mass is always a positive number.</p>
Modification	<p>Specifies the modifications from imported peptide files or from the Peptide Predictor tool.</p> <p>Available only for peptide databases.</p>
Num Isotopes Matched	<p>The number of isotopes matched in the expected calculated isotope spectra relative to the total number of isotopes used in the score calculation, in the format “x of y”, where</p> <ul style="list-style-type: none"> x = the number of isotopes matching the elemental composition used for the Isotopic Pattern Score calculation. y = the total number of isotopes considered in the Isotopic Pattern Score calculation. This is the number of isotope peaks expected to be above the spectral noise.
Peak Label	User-specified label.
Peptide Sequence	<p>Specifies the peptide sequence that the application uses to calculate the <i>m/z</i> or precursor <i>m/z</i>.</p> <p>Available only for peptide databases.</p>
PK	<p>Specifies if the peak was found. The application displays one of these flags:</p> <ul style="list-style-type: none"> A green circle when the peak is found A red square when the peak is not found <p>To display the peak results, hold your cursor over the indicator.</p> 
Protein Name	Alphanumeric name assigned to the protein.
Relative RT	Specifies the measured retention time relative to the RT Reference values set in the method.

Table 41. Common parameters for Compound Results and Sample Results tables (Sheet 8 of 8)

Column	Description
Response Ratio	The ratio of the Response value to the IS Response value. If the Response is specified as Area in the processing method, the units of both Response and IS Response are counts-sec. If the Response is specified as Height in the processing method, the units of both Response and IS Response are counts.
RSI/Rev Dot	Specifies the results of a reverse search of the NIST library. A reverse search compares a library entry to an unknown compound.
RT	Specifies the retention time for the compound.
RT Delta	Specifies the difference between the expected retention time and the measured retention time.
S/N	The signal-to-noise ratio calculated for the found peak.
Sample Amt	The injected volume multiplied by the conversion factor. For example, if you have 1000 ng/mL of a substance that is too concentrated for the mass spectrometer, you can dilute it by 1000. Then your injection volume is 1, your conversion factor is 1000, and your sample amount is 1000.
SI/Dot Product	Specifies the results of a forward search of the NIST library. A forward search compares the mass spectrum of an unknown compound to a mass spectral library entry.
Sample ID (Sample Results only)	A user-defined, alphanumeric string that identifies a sample.
Sample Name (Sample Results only)	A user-defined name that identifies a sample.
Sample Type (Sample Results only)	Defines how the application processes the sample data. Each sample is classified as one of the following sample types: Specimen, QC, Solvent, Calibrator, Hydrolysis, Unextracted, or Negative.
Std Add Amount	Specifies the standard amount added to the sample.
Theoretical Amt	Theoretical amount of the compound expected in the sample.
Type	In the Compound Results pane, specifies Target Compound, Confirming Peak, or Fragment. In the Sample Results pane, specifies Sample, Confirming Peak, or Fragment.

Inactive and Excluded Compounds

Use the Active and Excluded columns to control which compounds are used for calculating the calibration curve and for reporting.

Data Review - Batch_D [Quan]									
Compounds				Sample Results					
Flag	Compound	Comp		Flags	Flag Details		Active	Excluded	
▼	Aa	▼	Aa	▼	=	Aa	Aa	<input type="checkbox"/>	<input type="checkbox"/>
▶	FENTHION-CE2	Target Corr	▶	1			<input type="checkbox"/>	<input checked="" type="checkbox"/>	
▶	Sulfisomidine	Target Corr	▶	2			<input checked="" type="checkbox"/>	<input type="checkbox"/>	
				3			<input checked="" type="checkbox"/>	<input type="checkbox"/>	

Follow these procedures:

- [To make a sample active or inactive](#)
- [To exclude a calibration point](#)

❖ To make a sample active or inactive

1. Select the sample in the Sample Results pane.

All compounds in the selected sample appear in the Compounds pane. Inactive compounds are grayed out.

2. In the Compounds pane, select the compound whose active/inactive status you want to change.
 - When a compound is marked inactive, the application does not remove its data and calculated values from the result set. Instead, the application masks the appearance of that compound for that particular sample and grays the compound name in the compounds list.
 - When a calibration standard is marked inactive, the application no longer uses the data file's calibration point for the calibration and removes it from the graphical view of the calibration curve displayed in the Qualification pane. The calibration point is no longer part of the result set.

3. In the Sample Results pane, select or clear the **Active** check box.

Use the horizontal scroll bar at the bottom of the table to scroll to the Active column.

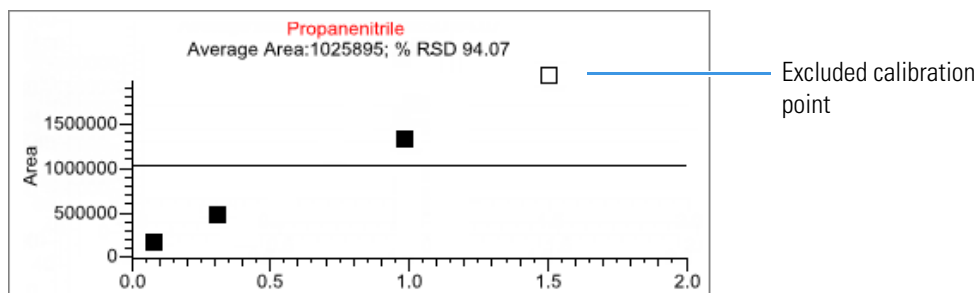
❖ To exclude a calibration point

In the sample list, select the **Excluded** check box for the sample.

Note Only calibration samples have the Excluded check box available. See [Inactive and Excluded Compounds](#).

Use the horizontal scroll bar at the bottom of the table to scroll to the Excluded column.

The application displays a value that is no longer used for calibration as an empty box in the graphical view of the calibration curve.



Exporting Compounds

You can export compound data to an Excel spreadsheet or to a CSV file. These export commands are available from the File menu in the Sample View, Compound View, and Comparative View.

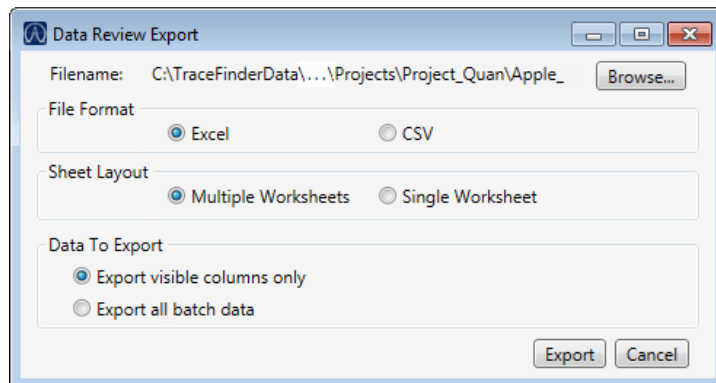
Follow these procedures:

- [To export compounds to an Excel spreadsheet](#)
- [To export compounds to a CSV file](#)

❖ To export compounds to an Excel spreadsheet

1. Choose **File > Export Data To > CSV or Excel** from the main menu.

The Data Review Export dialog box opens.



2. Click **Browse** and, in the Export Data to Excel dialog box, locate the folder where you want to save the file.

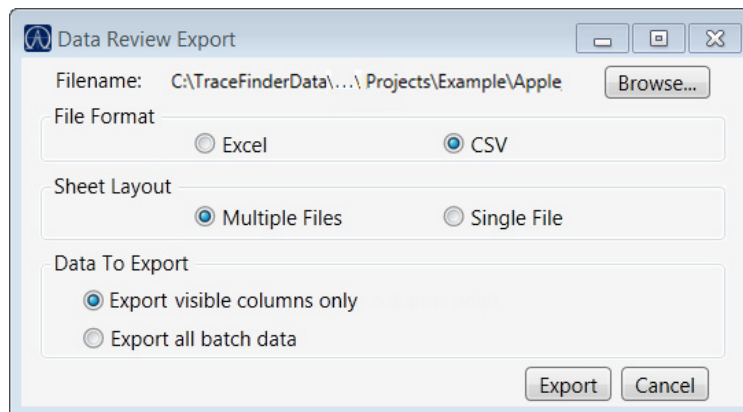
3. Type a file name for the XLSX file and click **Save**.
4. In the File Format area, in the Data Review Export dialog box, select the **Excel** option.
5. In the Sheet Layout area, select one of the following file formats for the spreadsheet.
 - **Multiple Worksheets:** Writes one sample to each Excel worksheet tab.
 - **Single Worksheet:** Writes all samples to a single Excel worksheet tab.
6. In the Data to Export area, select one of the following sets of data to export.
 - **Export Visible Columns Only:** Writes data from the displayed columns of selected samples to the specified worksheet format.
 - **Export All Batch Data:** Writes data from all columns (displayed or hidden) of all samples to the specified worksheet format.
7. Click **Export**.

The application saves the specified compound data to an Excel spreadsheet and opens the folder where you saved the file. The application names the file *Batch.xlsx*.

❖ To export compounds to a CSV file

1. Choose **File > Export Data To > CSV or Excel** from the main menu.

The Data Review Export dialog box opens.



2. Click **Browse** and, in the Export Data to Excel dialog box, locate the folder where you want to save the file.
3. Type a file name for the CSV file and click **Save**.
4. In the File Format area, in the Data Review Export dialog box, select the **CSV** option.
5. In the Sheet Layout area, select one of the following file formats for the spreadsheet.
 - **Multiple Files:** Writes one sample to each CSV file.
 - **Single File:** Writes all samples to a single CSV file.

6. In the Data to Export area, select one of the following sets of data to export.
 - **Export Visible Columns Only:** Writes data from the displayed columns of selected samples to the specified worksheet format.
 - **Export All Batch Data:** Writes all data from all samples in the batch to the specified worksheet format.
7. Click **Export**.

The application saves the specified compound data to a CSV spreadsheet.

When you selected to create multiple files, the application opens the folder where you saved the files. The application names each file *Batch_Compound.csv*.

When you selected to create a single file, the Excel application opens, displaying the exported data for all samples. The application names the file *Batch.csv*.

Working in the Report View for Quantitation Batches

The Report View displays example reports for the current batch. You must have an open batch to use the features in the Report View.

Follow these procedures:

- To open the Report View
- To preview a report
- To generate a report as a PDF, an Excel, or a CSV file
- To print a report
- To display a generated report
- To edit a report template
- To create a new report template

❖ To open the Report View

Click **Report View** in the navigation pane.



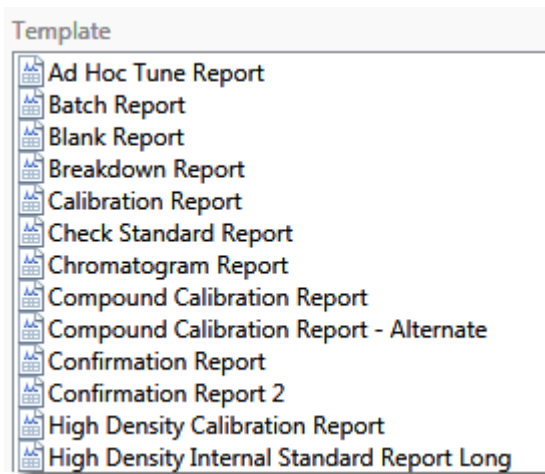
The application opens the [Report View](#).

❖ To preview a report

1. In the Template pane, select a report template.

The template list shows all the quantitation report templates that you configured in the Configuration console. Refer to “Specifying the Reports” in Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.

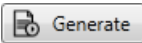
Figure 75. Example of a report template list



2. Click **Preview**, .

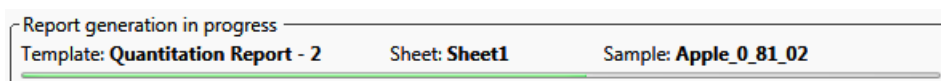
The application opens a Report Preview, showing the report information for the current batch in the selected report template format.

❖ To generate a report as a PDF, an Excel, or a CSV file

1. In the Template pane, select a report template.
2. Select the check box for each of the file types that you want to create: **PDF**, **Excel**, or **CSV**.
3. Click **Generate**, .

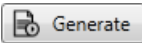
The application does the following:

- Displays a green progress bar as it generates the reports.



- Creates a report for the current batch as a PDF, an Excel, or a CSV file, using the selected report template format.
- Adds information about the generated report to the Generated Reports pane.
For details about the Generated Reports pane, see [Report View](#).
- Saves the report files to the ...\\TraceFinderData\\4.0\\Projects*batch*\\ReportOutput folder.

❖ To print a report

1. In the Template pane, select a report template.
2. Select the check box for the **Print** file format.
3. Click **Generate**, .

The application does the following:

- Creates a report for the current batch using the selected report template format.
- Prints the report to your default printer.
- Adds information about the generated report to the Generated Reports pane.

For details about the Generated Reports pane, see [Report View](#).

- Saves the report files to the ...\\TraceFinderData\\4.0\\Projects*batch*\\ReportOutput folder.

❖ To display a generated report

In the Generated Reports pane, click **View** for the report that you want to see.

Figure 76. Generated Reports pane showing a PDF report

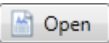
Generated Reports				
Template	Rule	Sample	Output	Generated Report File
Batch Report	Batch		pdf	View... Batch Report_20140113092939.pdf (3 pages)

Opens the generated file.

The application opens the output file.

❖ To edit a report template

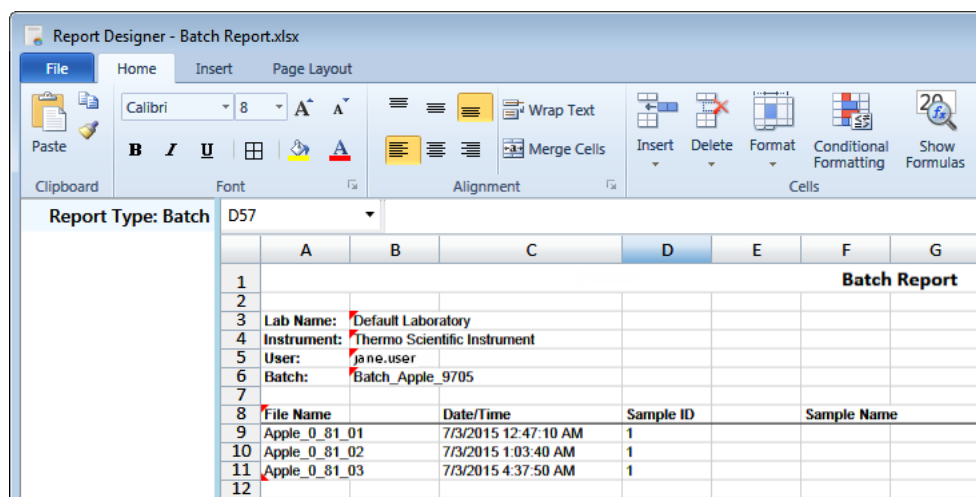
Note You cannot edit report templates that were provided with the TraceFinder application; however, you can open a TraceFinder template, make changes, and save it to a new template name.

1. In the Template pane, select a report template.
2. Click **Open**, .

The application opens the Report Designer showing the template in an Excel spreadsheet. See [Report Designer showing the template for the selected report](#).

Note When user security is activated, you must have Template Editing permission to edit report templates created by your laboratory. If the Open button is not active, user security is activated and you do not have Template Editing permission.

Figure 77. Report Designer showing the template for the selected report



Report Type: Batch		D57					
	A	B	C	D	E	F	G
1							
2							
3	Lab Name:	Default Laboratory					
4	Instrument:	Thermo Scientific Instrument					
5	User:	jane.user					
6	Batch:	Batch_Apple_9705					
7							
8	File Name	Date/Time	Sample ID				
9	Apple_0_81_01	7/3/2015 12:47:10 AM	1				
10	Apple_0_81_02	7/3/2015 1:03:40 AM	1				
11	Apple_0_81_03	7/3/2015 4:37:50 AM	1				
12							

3. Use the features in the Report Designer to edit the template.
See [Chapter 7, "Using the Report Designer."](#)
4. When you finish your changes, choose **File > Save** from the Report Designer menu bar.

❖ To create a new report template

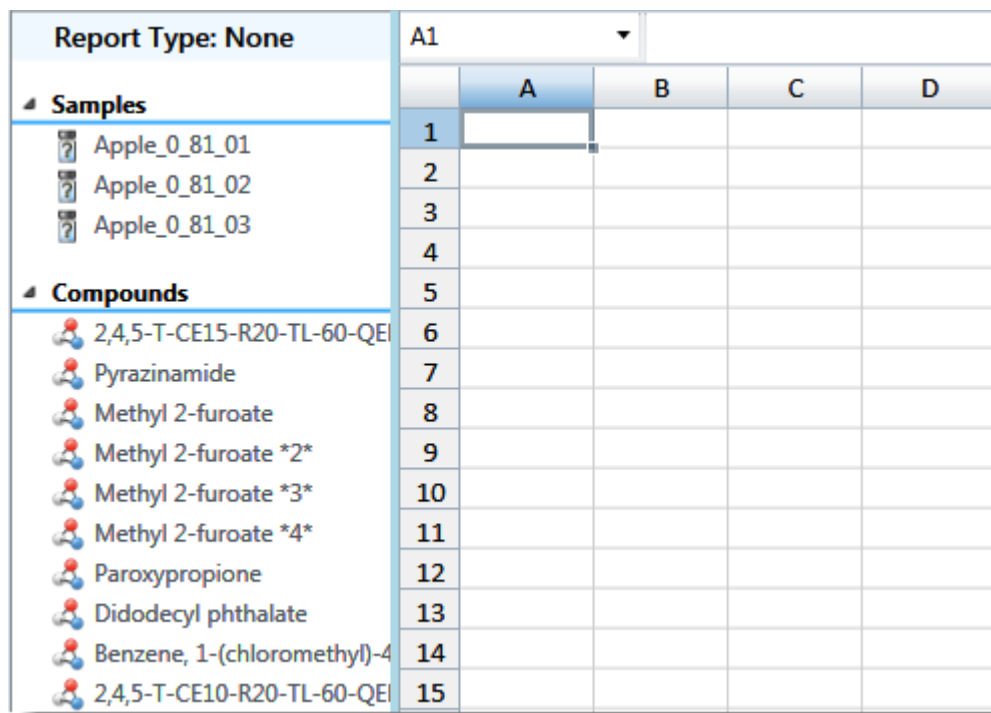
1. Click **New**, .

The application opens the [Report Designer](#) showing a new, empty template in an Excel spreadsheet.

The Report Type is None.

In the left pane, the spreadsheet lists all samples in the current batch and all compounds in the method used for the batch.

Figure 78. Report Designer showing a new, empty template



Report Type: None		A1			
		A	B	C	D
▲ Samples					
Apple_0_81_01					
Apple_0_81_02					
Apple_0_81_03					
▲ Compounds					
2,4,5-T-CE15-R20-TL-60-QEI					
Pyrazinamide					
Methyl 2-furoate					
Methyl 2-furoate *2*					
Methyl 2-furoate *3*					
Methyl 2-furoate *4*					
Paroxypropione					
Didodecyl phthalate					
Benzene, 1-(chloromethyl)-4					
2,4,5-T-CE10-R20-TL-60-QEI					

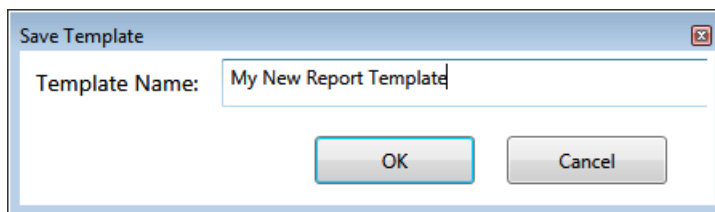
2. Use the features in the Report Designer to create the report template.

See [Chapter 7, “Using the Report Designer.”](#)

3. When you finish your changes, choose **File > Save** from the Report Designer menu bar.

The Save Template dialog box opens.

Figure 79. Save Template dialog box



Save Template

Template Name:

4. Type a name for the new report template and click **OK**.

Report View

Use the features in the Report View to display example reports for the current batch.

Figure 80. Report View

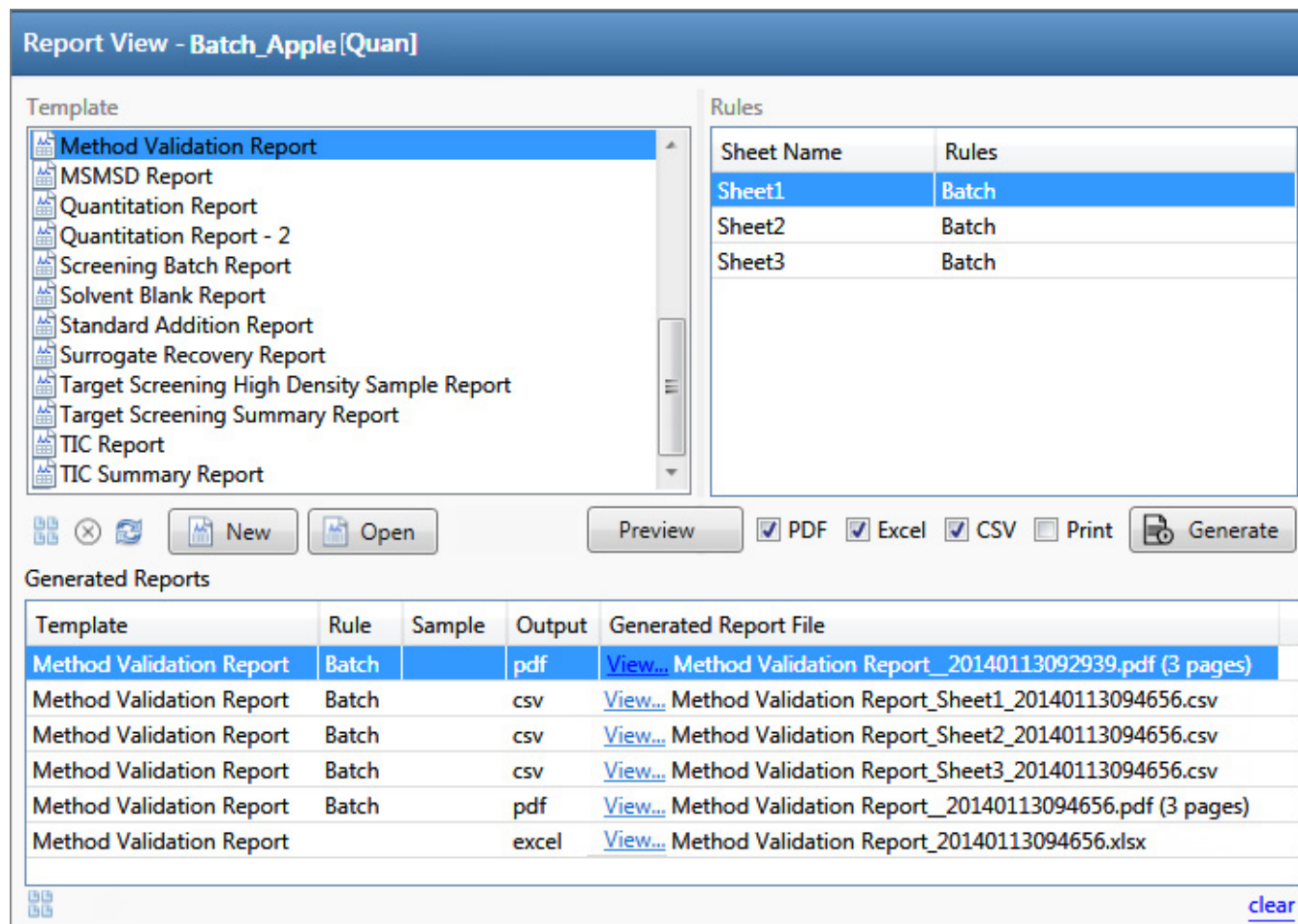

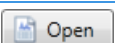


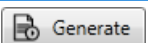



Table 42. Report View parameters (Sheet 1 of 2)

Parameter	Description
Template	
	Displays all report templates.
Rules	
Sheet Name	Specifies each sheet in the report.
Rules	Specifies the type of data used in each sheet in the selected report. <ul style="list-style-type: none"> • Batch • EachSample • SampleType: <i>SampleType</i> • CompoundType: <i>CompoundType</i> • SampleCustomFormula:

Table 42. Report View parameters (Sheet 2 of 2)

Parameter	Description
Buttons	
 View Report Templates	Displays the C:\TraceFinderData\4.0\Templates\ReportTemplates folder that contains all report templates.
 Open	Opens the selected report template in the Report Designer.
 New	Opens a blank report template in the Report Designer.
 Preview	Opens the Report Designer showing the report information for the current batch in the selected report template format.
PDF	Writes the generated report to a PDF file in the ...\\TraceFinderData\\4.0\\Projects\\ <i>batch</i> \\ReportOutput folder.
Excel	Writes the generated report to an Excel file in the ...\\TraceFinderData\\4.0\\Projects\\ <i>batch</i> \\ReportOutput folder.
CSV	Saves the generated report as a CSV file in the ...\\TraceFinderData\\4.0\\Projects\\ <i>batch</i> \\ReportOutput folder. When the report contains multiple sheets, the application writes each sheet as a separate CSV file.
Print	Prints the generated report to your default printer.
 Generate	Generates the selected type of reports for the current batch using the selected report template.
Generated Reports	
Template	Report template used for the report. See Example of a report template list .
Rule	Type of data used in each sheet of the report. See Rules .
Sample	For sample-level reports, the name of each sample in the report.
Output	Type of output specified for the report: PDF, Excel, CSV, or Print.
Generated Report File	Lists the output file name for each report in the ...\\TraceFinderData\\4.0\\Projects folder.
View	Displays the generated output file.
 View Generated Reports	Displays the C:\TraceFinderData\4.0\Projects folder that contains all report outputs.
Clear	Removes all reports from the Generated Reports display. This does not delete the reports from the C:\TraceFinderData\4.0\Projects folder.

Working in the Local Method View for Quantitation Batches

In the Local Method view, you can edit the local method parameters. A local method is a copy of a master method associated with a batch.

You can edit only the local copy of the method, or you can edit the master method and overwrite the local copy with the edited master method. Local methods are named *Batch_MasterMethod*.

❖ To open the Local Method View

1. Click **Analysis** in the navigation pane.
2. Click **Local Method**.



The Local Method view for the currently selected batch opens. See [Local Method view](#).

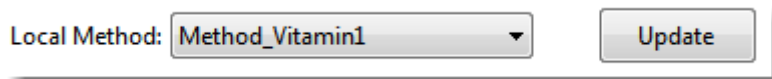
You can edit many of the method parameters in a local method. Editing the local method does not affect parameters in the master method.

For detailed descriptions of method parameters, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

3. Enter any local changes to the method.
4. When you have finished editing the local method, choose **File > Save**.
5. To process the batch or create new reports with the edited local method, return to the Batch View and submit the batch.

❖ To overwrite the local method with the master method in the Batch View

In the Batch View, click **Update**.



The application overwrites the local method with the master method of the same name. You can use this feature to overwrite an edited local method with the original master method or to overwrite the local method with an updated master method.

Figure 81. Local Method view

The screenshot displays the 'Local Method View' window for 'Equan1_408_Method_Equan_1*'. The left sidebar contains a navigation menu with the following items: 'Analysis' (expanded), 'Batch View', 'Data Review', 'Report View', 'Local Method' (selected), 'Acquisition', 'Quantitation', 'Processing', 'Compounds', 'QAQC', 'Groups', 'Intel Seq', 'Library Settings', and 'Reports'. The main panel shows the 'Master method: Method Equan_1' and various configuration fields: 'Lab Name' (Default Laboratory), 'Assay type' (Assay name), 'Injection Volume' (1.00), 'Mass Precision' (3.00), 'Ion range calc method' (Level), 'Use level' (dropdown), and 'Instrument Method' (Pesticides Using EQuan). 'Edit' and 'Update' buttons are located at the bottom right. A 'Notes' section with a text area is at the bottom left of the main panel.

Local Method View - Equan1_408_Method_Equan_1*	
Master method: Method Equan_1	
Lab Name:	Default Laboratory
Assay type:	Assay name
Injection Volume:	1.00
Mass Precision:	3.00
Ion range calc method:	Level
Use level:	
Instrument Method:	Pesticides Using EQuan
<div>Edit Update</div>	
Notes	

Using the Analysis Mode for Target Screening Batches

Use the features of the Analysis mode to do the following:

- Create target screening batches.
- Submit target screening batches for acquisition, processing, or report generation.
- Review target screening batches, batch data, reports, and local methods.

IMPORTANT TraceFinder 4.1 uses the same data as TraceFinder 4.0. By default, the application stores the method, project, template, and compound database data for the 4.1 release in the TraceFinderData\4.0 folder.

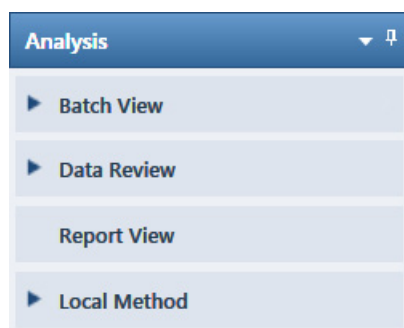
Contents

- [Working in the Batch View for Target Screening Batches](#)
- [Working in Data Review for Target Screening Batches](#)
- [Working in the Report View for Target Screening Batches](#)
- [Working in the Local Method View for Target Screening Batches](#)

❖ To access the Analysis mode

Click **Analysis** in the navigation pane.

The Analysis navigation pane opens.



Working in the Batch View for Target Screening Batches

In the Batch View, you can manually create and edit a new target screening batch or open and edit a previously saved batch. When you submit a batch, you can acquire and process data and optionally create reports for the submitted samples.

The Analysis mode includes a toolbar:

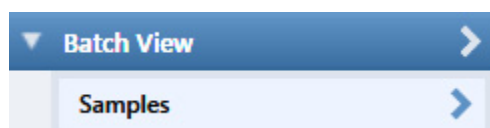


Use the [Toolbar](#) or the equivalent commands in the [Batch View Shortcut Menu](#) to create the sample list and submit samples for acquisition.

❖ To open the Batch View

1. Click **Analysis** in the navigation pane of the current mode.
2. Click **Batch View**.

The Batch View navigation pane opens.



The Batch View includes a single page of information, the Samples page.

Related Topics

- [Samples Page Features for Target Screening Batches](#)
- [Creating a New Target Screening Batch](#)
- [Editing a Target Screening Batch](#)
- [Submitting a Target Screening Batch](#)

Samples Page Features for Target Screening Batches

The Samples page is divided into two panes:

- Samples pane
Use the [Samples Pane](#) to create a batch.
- Compound Active Status pane
Use the [Compound Active Status Pane](#) to make specific compounds active or inactive.

Samples pane

Batch View - Batch_A [Screening]

Local Method: Method_Screening Update Instrument: Thermo Scientific Instrument User:

	Status	Filename	Groups	Blank Subtraction	Level	Sample ID	Sample name	Vial position	Injection volume	Conversion Factor	Channel	Barcode Expected	Barcode Actual	Sample Volume
▶ 1		Unknown1		<input type="checkbox"/>					10.0	1.000	Auto			1
2		Unknown2		<input type="checkbox"/>					10.0	1.000	Auto			1
3		Unknown3		<input type="checkbox"/>					10.0	1.000	Auto			1

▼▼▼▼

Compound Active Status

	RT	Compound	Active
▶ 1	1.91	FENTHION-CE20-R20-TL75-QED	<input checked="" type="checkbox"/>
2	2.72	Sulfisomidine	<input checked="" type="checkbox"/>

Compound Active Status pane

Tip To resize the panes, drag the separators that divide the panes.

Samples Pane

The samples pane includes the following features:

- [Column Display](#)
- [Status Indicators](#)
- [Groups](#)
- [Blank Subtraction](#)
- [Sample Weight Calculation](#)
- [Instrument Methods](#)
- [Toolbar](#)
- [Batch View Sample List](#)
- [Batch View Shortcut Menu](#)

Column Display

The sample list contains many columns of information. You can scroll to see all the columns, and you can customize which ones to display and their display order.

❖ To scroll the sample list

Use the horizontal scroll bar at the bottom of the sample list to view all the information.

When you use the scroll bar at the bottom of the sample list, the following columns remain fixed: Status, Filename, Groups, Blank Subtraction, Level, Sample ID, and Sample Name, while the other columns scroll right and left.

❖ To customize the column display

1. Right-click the sample list and choose **Modify Columns**.

The **Modify Columns dialog box** opens.

2. Use the arrow buttons to move all the columns that you want displayed to the Displayed Columns pane.

These columns appear after the Status, Filename, Groups, Blank Subtraction, Level, Sample ID, and Sample Name columns.

3. To arrange the order of the columns, do the following:

- a. In the Displayed Columns pane, select a column name.
- b. Use **Up** or **Down** to move the selected column up or down in the list.

The first column in the list represents the leftmost column in the Batch View sample list, and the last column in the list represents the rightmost column in the Batch View sample list.

Note The following columns are fixed: Status, Filename, Groups, Blank Subtraction, Level, Sample ID, and Sample Name.

4. To change the width of a column, do the following:

- a. In the Displayed Columns pane, select the column width.

5	Sample ID	100
▶ 6	Sample name	100
7	Vial position	100

- b. Type a new value for the width.

5. Repeat [step 4](#) for all columns whose widths you want to change, and click **OK**.

The columns in the sample list immediately reflect your changes. The application uses these settings for all sample lists in the Batch View.

Figure 82. Modify Columns dialog box

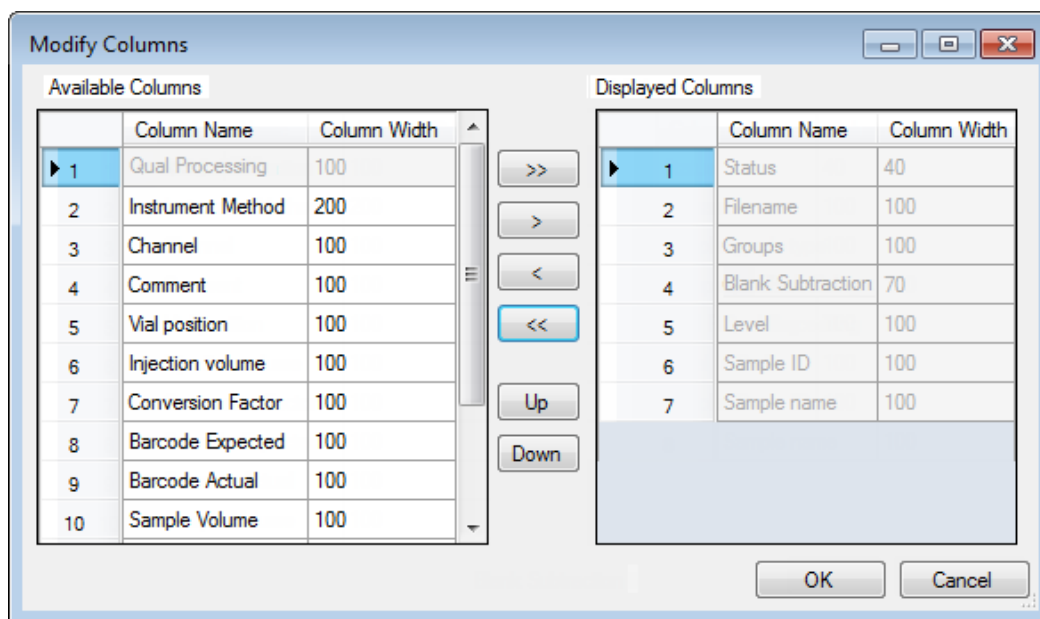




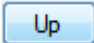
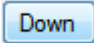






Table 43. Button descriptions for the Modify Columns dialog box




Button	Description
	Moves all columns to the Displayed Columns pane.
	Moves the selected column to the Displayed Columns pane.
The following buttons apply to all columns, except for those that are fixed: Status, Filename, Sample Type, Groups, Blank Subtraction, Level, Sample ID, and Sample Name.	
	Moves the selected column to the Available Columns pane.
	Moves all columns except fixed columns.
	Moves the selected column name in the Displayed Columns pane one row up in the column order.
	Moves the selected column name in the Displayed Columns pane one row down in the column order.

Status Indicators



Status indicators show the current status of each sample during the acquisition and processing.


-  Sample is not acquired.
-  Sample is acquired but not processed.
-  Sample is acquired and processed.
-  Sample is currently acquiring.

Local Method: Method_Screening Update

	Status	Filename	Groups	Level	Blank Subtraction	Sample ID	Sample name	V
1		Solvent_A			<input type="checkbox"/>			1
2		Cal_5			<input type="checkbox"/>			2
3		Unknown1			<input type="checkbox"/>			3

Status indicators

Note When you include unknown screening features in the target screening method and you choose to process with only the target screening criteria, the Samples view shows the Status for the samples as acquired and processed () , and the Unknown Screening View shows the Status for the samples as acquired but not processed () .

Note When you include unknown screening features in the target screening method and you choose to process with only the unknown screening criteria, both the Samples view and the Unknown Screening View show the Status for the samples as acquired and processed () .




Groups

Use the Groups feature to assign samples to a group.

❖ To create a group



- For each sample, type the name of a group in the Groups column.
Repeat this for each sample that you want to include in a group.
- Create as many groups as you want.

Note To assign a sample to multiple groups, separate the groups with a comma.

Status	Filename	Groups
	Benzo26473	groupB, groupA
	Benzo25557	groupB
	Benzo26154	groupB, groupA

Blank Subtraction

Use the Blank Subtraction feature to select which negative samples you want to use for peak subtraction. The application subtracts the areas of the peaks in the selected negative samples from the matching areas in the specimen samples.

Status	Filename	Groups	Level	Sample ID	Sample name	Blank Subtraction
	Apple_PosHCD_0_81_01			1		<input checked="" type="checkbox"/>
	Apple_PosHCD_0_81_02			1		<input type="checkbox"/>

When you process the batch sequence, the application subtracts the peaks in a selected negative sample from all specimen samples that follow it, until it encounters another negative sample.

To activate the Blank Subtraction feature, refer to “Editing the Processing Page” in Chapter 5, “Using the Method Development Mode for Target Screening Methods,” in the *TraceFinder Lab Director User Guide*.

Sample Weight Calculation

Use the sample weight features to calculate the conversion factor for a sample. The application uses different methods to calculate the conversion factor for liquid or solid calculation types.

Liquid: $SampleVolume \div DilutionFactor$

Solid: $(SampleVolume \times DilutionFactor) \div SampleWeight$

Manual: The application does not calculate the Conversion Factor. Instead, you can enter the Conversion Factor value.

Follow these procedures:

- To display the features for calculating sample weight
- To calculate the conversion factor for a liquid sample
- To calculate the conversion factor for a solid sample
- To manually specify the conversion factor for a sample

❖ To display the features for calculating sample weight

If the Conversion Factor, Sample Volume, Dilution Factor, Sample Weight, Calculation Type, and Final Units columns are not visible, right-click and choose **Enable Sample Weight Calculation**.

Conversion Factor	Sample Volume	Dilution Factor	Sample Weight	Calculation Type	Final Units
1.000	1	1	1	Liquid	
1.000	1	1	1	Solid	
1.000	1	1	1	Manual	

❖ To calculate the conversion factor for a liquid sample

Note The application uses the following formula to calculate the Conversion Factor:
 $SampleVolume \div DilutionFactor$

1. From the **Calculation Type** list, select **Liquid**.

For a liquid sample, the Sample Weight value is not editable.

2. In the Sample Volume column, type the volume in ng/mL for your sample.
3. In the Dilution Factor column, type the value for the dilution.

For example, if you have 1000 ng/mL of a substance that is too concentrated for the mass spectrometer, you can dilute it by 1000. Then your injection volume is 1, your conversion factor is 1000, and your sample amount is 1000.

4. In the Final Units column, type the units that you want to use for the calculated amount in the Data Review view or in reports.

❖ To calculate the conversion factor for a solid sample

Note The application uses the following formula to calculate the Conversion Factor:
 $(SampleVolume \times DilutionFactor) \div SampleWeight$

1. From the **Calculation Type** list, select **Solid**.
2. In the Sample Weight column, type the weight in ng for your sample.
3. In the Sample Volume column, type the volume in ng/ml for your sample.
4. In the Dilution Factor column, type the value for the dilution.

For example, if you have 1000 ng/ml of a substance that is too concentrated for the mass spectrometer, you can dilute it by 1000. Then your injection volume is 1, your conversion factor is 1000, and your sample amount is 1000.

5. In the Final Units column, type the units that you want to use for the calculated amount in the Data Review view or in reports.

❖ To manually specify the conversion factor for a sample

Note The application uses the specified conversion factor when it calculates the amount for the sample.

1. From the **Calculation Type** list, select **Manual**.

For a manually calculated sample, the only available columns are the Conversion Factor and the Final Units.


2. In the Conversion Factor column, type the conversion factor to use for your sample.
3. In the Final Units column, type the units that you want to use for the calculated amount in the Data Review view or in reports.

Instrument Methods

Use the Instrument Methods column to specify instrument methods for the samples.

Note By default, the Instrument Method column is not displayed in the Batch View sample list.




❖ To specify instrument methods for samples

1. Display the Instrument Method column in the sample list:
 - a. Right-click the sample list and choose **Modify Columns**.
The Modify Columns dialog box opens.
 - b. In the Available Columns pane, select **Instrument Method**.
 - c. Click  to move the Instrument Method column to the Displayed Columns pane.
 - d. Click **OK**.

The application displays the Instrument Method column, defaulting to the instrument method specified in the master method.

2. Click the Instrument Method column and select an instrument method from the list.
This list contains all the available instrument methods. Instrument methods from external sources are prefixed with “Ext:”.

You can specify a different instrument method for each sample.

Batch View - Batch_8266 [Screening]						
Local Method: Method_Screening		Update		Instrument: Thermo Scientific Instrument		
	Status	Groups	Filename	Sample type	Blank Subtraction	Instrument Method
4			Unknown4	Specimen	<input type="checkbox"/>	Ext: Instrument1
5			Unknown5	Specimen	<input type="checkbox"/>	Ext: Instrument2
6			Unknown6	Specimen	<input type="checkbox"/>	Ext: Instrument3

When you submit the batch for acquisition, the application saves a copy of the selected instrument methods to the following folders:

External instrument methods:

...\\TraceFinderData\\4.0\\Projects\\...\\batch\\Methods\\method\\ExternalMethods

Local instrument methods:

...\\TraceFinderData\\4.0\\Projects\\...\\batch\\Methods\\method

Toolbar

The Analysis mode includes this toolbar for creating and submitting a batch.



Table 44. Toolbar icons

Icon	Description
	Adds the specified number of new, empty samples to the end of the sample list. See the instructions To add samples to the list .
	Inserts a new, empty sample or samples above the selected sample. See the instructions To insert samples into the list .
	Removes the selected samples from the sample list. See the instructions To remove samples from the list .
	Adds imported samples from a CSV, an XML, or an SLD file to the sample list. See the instructions To import samples into the list .
	Submits only the selected samples for acquisition, processing, or report generation. See the instructions To submit samples in the batch .
	Submits the batch for acquisition, processing, or report generation. See the instructions To submit samples in the batch .
	Submits only the selected samples for processing. See the instructions To submit selected peaks for processing .
	Opens the Acquisition mode where you can use a batch template to define a standard sequence composed of various sample types to be assembled into a batch of samples. See Working in Data Review for Target Screening Batches .
	Opens the Acquisition mode where you can create a batch template that contains the basic settings and sample types for your batches. See Using the Acquisition Mode .
	Opens the Quick Acquisition window where you can quickly submit a single sample. See Appendix A, "Using Quick Acquisition."
	Opens the Audit Viewer where you can view audit logs. See Chapter 8, "Using the Audit Viewer." Available only when you enable Auditing in the Administrator Console. Refer to the instructions in the <i>TraceFinder Administrator Console User Guide</i> .

Batch View Sample List

The sample list displays all the quantitative data for the samples of a batch.

Status indicators for each sample indicate if the sample is currently acquiring, not acquired, acquired, or processed.

The sample list includes the following columns of information:

Figure 83. Batch View sample list




Local Method:

Method_Screening

Update

Instrument:

Thermo Scientific Instrument

	Status	Filename	Groups	Level	Sample ID	Sample name	Blank Subtraction	Vial position	Injection volume
1		UnknownA1			1		<input type="checkbox"/>	CStk1-01:7	10.0
2		UnknownA2			1		<input type="checkbox"/>	CStk1-01:7	10.0
3		UnknownA3			1		<input type="checkbox"/>	CStk1-01:7	10.0

Calculation Type	Conversion Factor	Dilution Factor	Sample Weight	Sample Volume	Final Units	Sample type
Liquid	1.000	1	1	1		Specimen
Liquid	1.000	1	1	1		Specimen
Liquid	1.000	1	1	1		Specimen

Instrument Method	Channel	Barcode Expected	Barcode Actual	Comment
Instrument1	Auto			
Instrument1	Auto			
Instrument1	Auto			

Table 45. Batch View sample list columns (Sheet 1 of 3)





Column	Description
Status	 Sample is not acquired.  Sample is acquired but not processed.  Sample is acquired and processed.  Sample is currently acquiring.
Filename	Name of the raw data file that contains the sample data.
Groups	Threshold group to which a sample belongs.
Blank Subtraction	Specifies a negative sample to use for blank subtraction.
Level	The level defined for a calibration sample or quality control sample.

Table 45. Batch View sample list columns (Sheet 2 of 3)

Column	Description
Sample ID	A user-defined, alphanumeric string that identifies a sample.
Sample Name	A user-defined name that identifies a sample.
Vial Position	The tray vial number used for an autosampler acquisition.
Injection Volume	<p>The injection volume (in microliters) of the injected sample.</p> <p>When you are using an autosampler, you can set the default injection volume in the Autosampler dialog box in the Instrument View. The minimum and maximum injection volumes that you can use depend on the Autosampler you configure. The usable range depends on the injection mode and might be smaller than the displayed range.</p> <p>The Injection Volume value set in the master method overwrites the value in the instrument method.</p> <p>Valid range: 0.1 through 5000 µL</p>
Calculation Type	<p>Liquid: The application calculates the Conversion Factor as</p> $\text{SampleVolume} \div \text{DilutionFactor}$ <p>Solid: The application calculates the Conversion Factor as</p> $(\text{SampleVolume} \times \text{DilutionFactor}) \div \text{SampleWeight}$ <p>Manual: Sample Volume, Dilution Factor, Sample Weight, and Final Units columns are not available, and the Conversion Factor value is editable.</p>
Conversion Factor	<p>Editable only when Calculation Type is Manual.</p> <p>Default: 1</p>
Sample Volume	Default: 1
Dilution Factor	Default: 1
Sample Weight	<p>Available only when Calculation Type is Solid.</p> <p>Default: 1</p>
Final Units	<p>Specifies the calculated amount in the Data Review view or in reports.</p> <p>Default: 1</p>
Instrument Method	Specifies the instrument to use for the acquisition. This column is hidden by default. To display this column, see To customize the column display .
Channel	Specifies the channel on which the sample was run. If the sample is not acquired, the value is Pending. The Channel column is available only when you have activated multiplexing in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the <i>TraceFinder Lab Director User Guide</i> .

Table 45. Batch View sample list columns (Sheet 3 of 3)

Column	Description
Barcode Expected	A user-entered barcode for the vial.
Barcode Actual	An actual barcode for the vial. This value is not editable.
Comment	A user-defined comment for the sample.
Sample Type	Defines how the application processes the sample data. Each sample is classified as one of the following sample types: Specimen, QC, Solvent, Calibrator, Hydrolysis, Unextracted, or Negative. Default: Specimen


Batch View Shortcut Menu

The Batch View includes a shortcut menu for creating a batch.

Table 46. Batch View shortcut menu commands (Sheet 1 of 2)

Command	Description
Add Sample	Adds a single empty row to the sample grid.
Insert Sample	Inserts a single empty row to the sample grid above the selected row.
Insert Copy Sample	Copies the currently selected row and inserts a copy above the row.
Reinject Selected Samples	Creates a copy of the selected sample and appends INJ001 to the file name. Additional reinjections of the same sample are numbered INJ002, INJ003, and so forth.
Remove Selected Samples	Removes selected samples from the sample grid.
Import Samples	Opens the Sample Import Tool. See To import samples into the list .
Browse in Raw File (Move)	Opens a dialog box where you can select a raw data file to use for the selected sample row. The application removes the raw data file from the source location.
Browse in Raw File (Copy)	Opens a dialog box where you can select a raw data file to use for the selected sample row. The application copies the raw data file from the source location.
Map Raw Files to Samples	Opens a dialog box where you can select multiple raw data files to use for the selected sample rows.
Copy Down	Copies the value in the selected row to all rows below it. This command is available only when you have selected a value that can be copied down.
Fill Down	Enters sequential values in the column starting with the value in the selected row and ending with the last row in the column. This command is available only when you have selected a value that can be filled down.
Modify Columns	Opens the Modify Columns dialog box .
Hide Sample Weight Calculation	Displays or hides the Sample Volume, Dilution Factor, Sample Weight, Calculation Type, and Final Units columns.

Table 46. Batch View shortcut menu commands (Sheet 2 of 2)

Command	Description
Copy	Copies the data in the selected rows or columns to the Clipboard. Use this command to copy sample information into a text editor or spreadsheet application. You cannot paste this data back into the Batch View sample list.
Copy with Headers	<p>Copies the data in the selected rows or columns and the associated column headers to the Clipboard. Use this command to copy sample information into another text editor or spreadsheet application. You cannot paste this data back into the sample list.</p> <p>For example</p>  <p>Copy With Headers from TraceFinder</p> <p>Paste into Excel spreadsheet</p>
Paste	Pastes a single column of copied data from another text editor or spreadsheet application into the selected column.
Undo Last Paste	Removes the last pasted item in the Batch View.
Export to CSV File	Opens the Save As dialog box where you can save the current sample list to a CSV file.
Edit Instrument Method	<p>Opens the Instrument Setup window where you can edit the parameters of the instrument method.</p> <ul style="list-style-type: none"> • When you edit an external method, the application updates the method in the ...\\Xcalibur\\methods folder. • When you edit an internal method, the application updates the method in the ...\\TraceFinderData\\4.0\\Projects\\...\\batch\\Methods\\method folder. <p>For detailed information about editing instrument methods, refer to Chapter 3, “Using Instrument Methods in the Method Development Mode,” in the <i>TraceFinder Lab Director User Guide</i>.</p>
View Sample in Qual Explorer	Displays the selected sample in the qualitative explorer application that you configured as your default in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the <i>TraceFinder Lab Director User Guide</i> .

Compound Active Status Pane

In the Compound Active Status pane, you can choose specific compounds to be active or inactive.

❖ To set a compound as active or inactive

1. In the sample list, select a sample.

All compounds in the selected sample are listed in the Compound Active Status pane.

Compound Active Status

	RT	Compound	Active
▶ 1	1.91	FENTHION-CE20-R20-TL75-QED	<input checked="" type="checkbox"/>
2	2.72	Sulfisomidine	<input checked="" type="checkbox"/>

The default active/inactive status is determined by the identification settings in the local method. For information about setting the identification parameters, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

- To display compounds alphabetically, right-click and choose **Sort by Compound Name**.
 - To display compounds from shorter to longer retention times, right-click and choose **Sort by Retention Time**.
2. Select or clear the **Active** check box for the compound.

Compound Active/Inactive Status

You can specify which compounds are active or inactive in the [Local Method View](#) or the [Batch View](#).

Figure 84. Active and inactive compounds in the Local Method View



Local Method View - Batch_1_Method_Benzos					
Master method: Method Benzos					
Acquisition List	Identification	Detection	Calibration	Calibration levels	QC Check levels
	RT	Compound	Compound type	Active	CAS No
1	1.91	FENTHION-CE20-R20-TL...	Target Compound	<input type="checkbox"/>	55389
▶ 2	2.72	Sulfisomidine	Target Compound	<input checked="" type="checkbox"/>	515640

For details about setting the status on the Identification page, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

5 Using the Analysis Mode for Target Screening Batches

Working in the Batch View for Target Screening Batches

Figure 85. Active and inactive compounds in the Batch View

Batch View - Batch_2 [Screening]						
Local Method: Method_Screening				Update		
	Status	Filename	Groups	Level	Blank Subtraction	Sample ID
▶ 1		VitWaterEquanMaxA001			<input type="checkbox"/>	CStk1-01:3
2		VitWaterEquanMaxA002			<input type="checkbox"/>	CStk1-01:4
Compound Active Status						
	RT	Compound	Active			
▶ 1	5.88	1-Butanol, 4-(butylnitrosoami...	<input type="checkbox"/>			
2	16.68	3,3'-Diaminodiphenylmethane	<input checked="" type="checkbox"/>			
3	2.91	4-Methyl-2,6,7-trioxa-1-phosp...	<input checked="" type="checkbox"/>			

For details about setting the status in the Batch View, see [Compound Active Status Pane](#).

Creating a New Target Screening Batch

In the Batch View, you can create a new batch.

Follow these procedures:

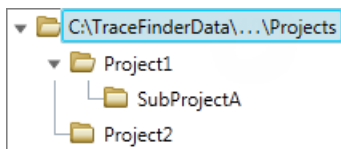
- [To create a new batch](#)
- [To add samples to the list](#)
- [To insert samples into the list](#)
- [To import samples into the list](#)
- [To remove samples from the list](#)
- [To copy a sample](#)
- [To reinject a sample](#)
- [To edit sample values](#)
- [To browse in raw data files](#)

❖ To create a new batch

1. Choose **File > New > Batch** from the main menu.

The [Create New Batch Dialog Box](#) opens and displays all drives that contain projects.

2. Select a drive from the list.



Tip The application displays all configured and enabled repositories.

3. Select the folder where you want to store your batch.

Tip To activate the Create button, you must enter a unique batch name. If the Create button is not activated, you have entered a batch name that is already used.

To create a new folder for the storage location, see [Editing Folders for Batches](#).

4. Select **Screening** from the Type list.

The batch list displays all batches in the selected folder. The Method list displays all methods for the target screening type.

5. Select a master method from the Master Method list.

5 Using the Analysis Mode for Target Screening Batches

Working in the Batch View for Target Screening Batches

6. Click **Create**.

A new batch opens with one Specimen sample.

Batch View - Batch_13 [Screening]						
Local Method: Method_Screening					Update	
	Status	Filename	Groups	Level	Blank Subtraction	Sample ID
▶1		Unknown1			<input type="checkbox"/>	CStk1-01:3

The batch name in the title bar indicates that you are creating a target screening batch.

❖ To add samples to the list

1. To add a single sample row, right-click the sample list and choose **Add Sample**.
2. To add multiple sample rows, select the number of rows and then click the **Add Sample** icon, .

The application adds the specified number of new, empty samples to the end of the sample list.

❖ To insert samples into the list

Select the sample above which you will insert new, Specimen samples, and then do one of the following:

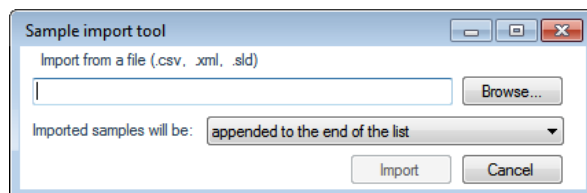
- To insert a single sample row, right-click and choose **Insert Sample**.
- To insert multiple sample rows, select the number of rows and then click the **Insert Sample** icon .

The application inserts the Specimen samples above the selected sample.

❖ To import samples into the list

1. Choose **Batch > Import Samples** from the main menu, or click the **Import Samples** icon, .

The Sample Import Tool dialog box opens.



From this dialog box, you can import samples from a CSV, an XML, or an SLD file.

2. Click **Browse** and select a CSV, an XML, or an SLD file that contains the samples to import.

- From the Imported Samples Will Be list, select **Appended to the End of the List** or **Inserted at the Selected Row**.
- Click **Import**.

The Sample Import Tool dialog box closes, and the application adds the specified samples to the sample list.

When you import samples from an Xcalibur sequence file (.sld), the application makes the following column name substitutions:

Xcalibur column	TraceFinder column
Position	Vial Position
Inj Vol	Injection Volume
Dil Factor	Conversion Factor

When you import samples from an Xcalibur sequence file (.sld), the application makes the following sample type substitutions:

Xcalibur sample type	TraceFinder sample type
Blank	Negative
Std Bracket	Calibrator

- (Optional) When using multiplexing, select a channel for each imported sample.
Imported samples default to Auto.

Note The Channel column is available only when you have activated multiplexing in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.

❖ To remove samples from the list

- Select the samples that you want to remove.

Tip Use the CTRL or SHIFT keys to select multiple samples.

- Right-click and choose **Remove Selected Samples**.

❖ To copy a sample

- Select the sample that you want to copy.
- Right-click and choose **Insert Copy Sample**.

The application inserts the copy above the selected sample.

❖ To reinject a sample

1. In the sample list, select the sample that you want to reinject.
2. Right-click and choose **Reinject This Sample**.

The application creates a copy of the selected sample and appends INJ001 to the file name. Additional reinjections of the sample are numbered INJ002, INJ003, and so forth. The application copies all parameter values from the original sample.

❖ To edit sample values

1. For each sample, do one of the following:

Type a new file name over the current filename.

—or—

Double-click the Filename column and locate a raw data file to use for the sample.

—or—

Right-click and choose **Browse in Raw File**, and then locate a raw data file to use for the sample.

By default, the application sets the Sample Type to Unknown.

2. For each sample, click the Sample Type column and select a sample type from the list.

Available sample types

Specimen	Hydrolysis	Solvent	QC
Unextracted	Calibrator	Negative	

3. For each Calibrator or QC sample, select a level from the Level list.

The sample levels are defined in the master method. If there are no levels to select in the Level list, do the following:

- a. Return to the Method Development mode.
- b. Open the method.
- c. Click the **Compounds** tab.
- d. Click the **Calibration Levels** tab.
- e. Add the levels.
- f. Save the method.
- g. Return to the Batch View in the Analysis mode, and then click **Update**.

Local Method: Method_Alprazolam_Screening ▼

Update

The application updates the local method with the new sample levels.

For detailed instructions about specifying calibration levels, refer to Chapter 4, “Using the Method Development Mode for Quantitation Methods,” in the *TraceFinder Lab Director User Guide*.

4. Type a vial position in the Vial Position column for each sample.
5. Type a volume in the Injection Volume column for each sample.

The minimum injection volume value allowed is 0.1 µL; the maximum injection volume value allowed is 5000 µL.

6. (Optional) Type or edit the values for the remaining columns.

Note When you use the scroll bar at the bottom of the sample list, the following columns remain fixed: Status, Filename, Sample Type, Groups, Blank Subtraction, Level, Sample ID, and Sample Name, while the other columns scroll right and left.

To automatically copy or fill values in these columns, see [Appendix B, “Using Copy Down and Fill Down.”](#)

❖ To browse in raw data files

1. Do one of the following:

Double-click the Filename column.


—or—

Right-click and choose **Browse in Raw File**.

The What Raw File Would You Like to Use dialog box opens.

2. Select a raw data file to use for the sample or use the CTRL key to select multiple files, and then click **Open**.

The application overwrites the selected, unacquired sample in the batch with the first “browsed in” file and adds any additional browsed in files below the selected sample.

For all browsed-in raw data files, the application sets the Status to Acquired, , and sets the Sample Type to Specimen.

Note You cannot overwrite an acquired sample. When you select a sample that is acquired, the application adds all browsed in files below the selected sample.

5 Using the Analysis Mode for Target Screening Batches

Working in the Batch View for Target Screening Batches

Create New Batch Dialog Box

Use the Create New Batch dialog box to select a folder and method for your batch and to name the new batch.

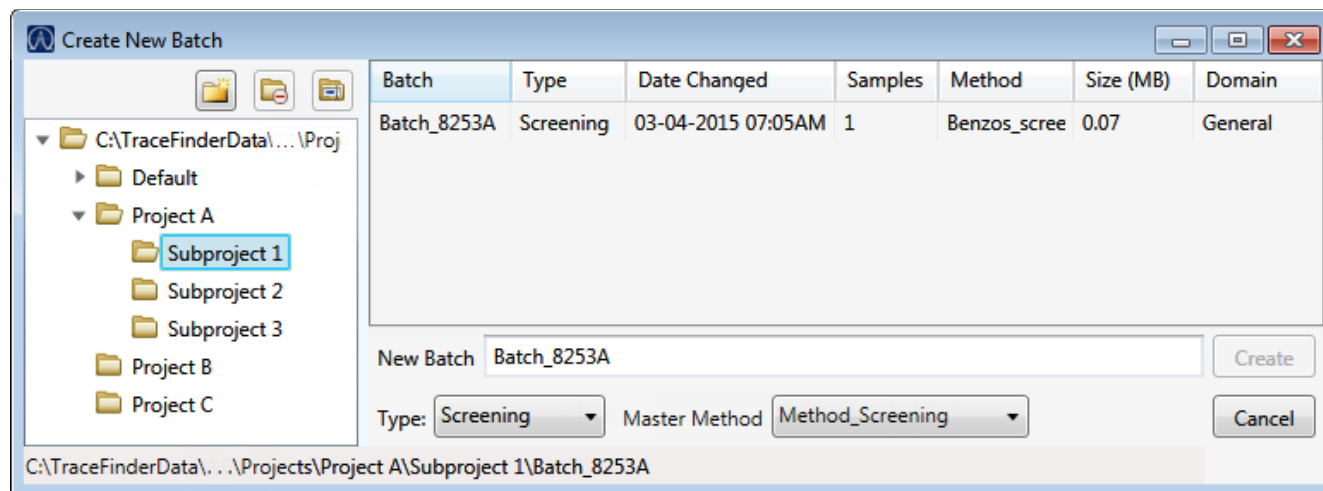


Table 47. Create New Batch dialog box parameters (Sheet 1 of 2)




Parameter	Description
 Create New Folder	Adds one of the following: <ul style="list-style-type: none">• When a drive is selected, adds a new project-level folder to the drive.• When a project folder is selected, adds a subproject-level folder to the selected project.• When a subproject folder is selected, adds a lower-level folder to the subproject. Or, you can right-click and choose Create Folder.
 Delete Folder	With no confirmation prompt, immediately removes the selected folder. You cannot delete a folder that contains lower-level folders; you must delete the lower-level folders first. Or, you can right-click and choose Delete.
 Rename Folder	Renames the selected folder. Or, you can right-click and choose Rename.
Batch table	
Batch	Name of batches in the selected project.
Type	Type of batch: Quan, Screening, or Unknown Only.
Date Changed	Date that the batch was last updated.
Samples	Number of samples in the batch.
Method	Name of the method used to create the batch.
Size	Size of the batch in megabytes.

Table 47. Create New Batch dialog box parameters (Sheet 2 of 2)

Parameter	Description
Domain	TraceFinder domain in which the batch was created.
New batch parameters	
New Batch	Name of the new batch to create. Note If the Create button is not activated, you have entered a name that is already used or you have not selected a method.
Type	Type of batch to create: Quan, Screening, or Unknown Only.
Method	Method used to create the new batch.
Path	Path to the project in the TraceFinderData\4.0\Projects folder where the batch is created.
Buttons	
Create	Creates the specified batch and opens the Batch View for the new batch.
Cancel	Closes the Create New Batch dialog box without creating a batch.


Editing Folders for Batches

From the Create New Batch dialog box, you can create new folders for your batches. You can also delete or rename folders.

Use these procedures:

- [To create new project folders](#)
- [To delete project folders](#)
- [To rename project folders](#)

❖ To create new project folders


1. In the Create New Batch dialog box, select the folder for which you will create a new lower-level folder.
 - You can select the main TraceFinderData\4.0\Projects folder and create a new folder under it.
 - You can select one of the existing folders and create a lower-level folder under it.
2. Click the **Create Folder** icon, .

The application adds a new lower-level folder to the selected folder.
3. Select the new folder name and type a name for the folder.

Folder names are limited to 30 characters and can contain spaces and special characters, except for the following special characters: \ / : + ? " < >

Note After you add a lower-level folder, you cannot rename the parent folder.


❖ To delete project folders

1. In the Create New Batch dialog box, select the folder to delete.
2. Click the **Delete Folder** icon, .

With no confirmation prompt, the application immediately removes the selected folder.

Note You cannot delete folders that contains lower-level folders; you must delete the lower-level folders first.

❖ To rename project folders

1. In the Create New Batch dialog box, select the folder to rename.
2. Click the **Rename Folder** icon, .

Note You cannot rename folders that contain lower-level folders.

3. Type a new name for the folder and press ENTER.

The application saves the new folder name.

Editing a Target Screening Batch

In the Batch View, you can open a saved batch and edit the sample list. You can add samples, edit samples, or remove samples. If the batch has already been acquired, you can select specific samples for reinjection. If the batch has unacquired samples when you complete your edits, you can save it as a “ready to acquire” batch.

Follow these procedures:

- [To open a saved batch](#)
- [To open a recent batch](#)
- [To edit samples in a batch](#)
- [To reinject a sample from a previously acquired batch](#)

❖ To open a saved batch

1. Choose **File > Open > Batch** from the main menu.

The [Open Batch dialog box](#) opens.

2. Select a project and a subproject.
3. Select **Screening** or **Any** from the Type list.

The batch list displays all batches created with Target Screening methods (or when you select Any, all methods of all types).

4. Select a batch from the list.

5. Click **Open**.

The selected batch opens in the Batch View.

Figure 86. Open Batch dialog box

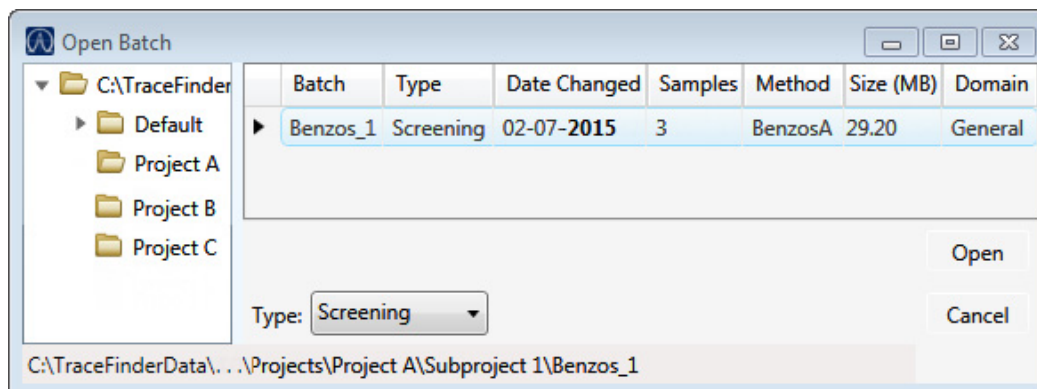


Table 48. Open Batch dialog box parameters

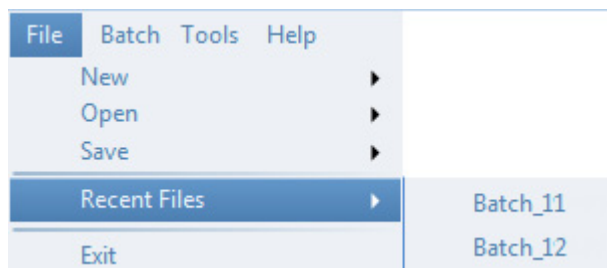
Parameter	Description
Batch	Name of batches in the selected project.
Type	Type of batch: Quan, Screening, or Unknown Only.
Date Changed	Date the batch was last updated.
Samples	Number of samples in the batch.
Method	Name of the method used to create the batch.
Size	Size of the batch in megabytes.
Domain	TraceFinder domain in which the batch was created.
<i>Path</i>	Path to the project in the TraceFinderData\4.0\Projects folder where the batch is stored.
Buttons	
Type	Type of batch to display in the Batch list: Quan, Screening, Unknown Only, or Any.
Open	Opens the Batch View for the selected batch.
Cancel	Closes the Open Batch dialog box without opening a batch.

5 Using the Analysis Mode for Target Screening Batches

Working in the Batch View for Target Screening Batches

❖ To open a recent batch

Choose **File > Recent Files > *batch*** from the main menu.



The selected batch opens in the Batch View.

❖ To edit samples in a batch

Use the commands described in [Working in the Batch View for Target Screening Batches](#).





You can add new samples, edit samples, or delete samples.

❖ To reinject a sample from a previously acquired batch

1. In the sample list, select the sample that you want to reinject.
2. Right-click and choose **Reinject This Sample**.

The application creates a copy of the selected sample and appends INJ001 to the file name. Additional reinjections of the sample are numbered INJ002, INJ003, and so forth. The application copies all parameter values from the original sample.

A green status icon indicates previously acquired samples (acquired and processed), and the sample name is grayed out. A blue status icon indicates samples created for reinjection (not acquired).

	cal_50_INJ001	Calibrator	10
	cal_50	Calibrator	10
	cal_10_INJ001	Calibrator	10
	cal_10	Calibrator	10

When you submit all samples in this batch, the application acquires all samples (including previously acquired samples).

3. To save this batch with the new samples for reinjection, choose **File > Save > Batch** from the main menu.

The batch is saved as a prepared batch that is ready to submit. You can open this batch from the Reinject Samples page in the Acquisition mode and submit the batch. The application acquires only the samples that have not been previously acquired.

Submitting a Target Screening Batch


In the Batch View, you can submit an entire batch or only selected samples in the batch. When you submit a batch for acquisition and processing, you can choose to create reports for the submitted samples. See [Submit Options dialog box](#).

For a description of commands in the shortcut menu, see [Batch View shortcut menu commands](#).

Follow these procedures:



- [To submit selected peaks for processing](#)
- [To submit samples in the batch](#)
- [To view the output files](#)

❖ To submit selected peaks for processing

1. In the Peak List, select the **Selected** check box for the peaks that you want to process.
2. Click the **Submit ... for Processing** icon, .

The application processes the selected peaks and updates the data in the Peak Identifications pane.

❖ To submit samples in the batch

1. Do one of the following:
 - To submit all samples in the batch, click the **Submit Batch** icon, .
 - To submit specific samples, select the samples and click the **Submit Selected Samples** icon, .

The [Submit Options dialog box](#) opens.

Note You can also submit only selected peaks for processing. See [To submit selected peaks for processing](#).

2. To acquire (or reacquire) the submitted samples, select the **Acquire Data** check box.
 - When all submitted samples have been previously acquired, this option is (by default) not selected.
 - When one or more samples in the batch have not been acquired, this option is (by default) selected.

Tip You can process the data with or without performing peak detection. For example, you might want to turn off peak detection when reprocessing samples.

3. To process the submitted samples, select the **Process Data** check box.

The application displays processing options for the Target Screening method.

If the method includes unknown screening features, the application also displays unknown screening options. For specifying processing data for batches that also use unknown screening features, see [To submit samples in the batch](#) in Chapter 6, “Using the Analysis Mode for Unknown Screening Batches.”

4. Select the check box for the options that you want to use.

Peak Detect: Performs peak detection. You can process the data with or without performing peak detection. For example, you might want to turn off peak detection when reprocessing samples.

Identify and Confirm: Performs both identification and confirmation.

5. (Optional) Select the **Create Reports** check box.
6. (Optional with multiplexing activated) Select the **Priority Sequence** check box.

The application acquires the priority batch on the next available channel or the assigned channel.

7. (Optional without multiplexing activated) Select the **Priority Sequence** check box and then select one of the following priority options to place the batch in the queue:
 - **Next Available Batch** places the batch immediately after the currently acquiring batch.
 - **Next Available Sample** places the batch immediately after the currently acquiring sample.

Note When you select Full Sequence Submission in the Configuration console, these options are unavailable because the current batch and the current sample are, in effect, the same thing.

8. To specify the following optional parameters, click **Show Details**.
 - a. Select the **Use** check box for the device that you want to use for this acquisition.
 - b. Select the **Start Device** check box to indicate the device that will initiate communication with the other instruments.

This is usually the autosampler.
 - c. Select the **Start When Ready** check box, which starts all instruments together when they are all ready.

When this is cleared, individual instruments can start at different times and then have to wait for the last instrument to be ready.
 - d. Select the system state after it acquires the last batch: **On**, **Standby**, or **Off**.
9. To start the selected processes, click **OK**.

The selected processes begin, and the application shows the real-time display at the bottom of the current window. You can begin another batch in the Analysis mode while you watch the real-time display of the currently acquiring batch.

IMPORTANT When your batch also uses unknown screening features, after the first processing, you might see an error message stating that the total number of results is too large for a single batch (more than 500 000). Return to the Processing pages for the unknown screening features (refer to Chapter 5, “Using the Method Development Mode for Unknown Screening Methods,” in the *TraceFinder Lab Director User Guide*) and make one or more of these parameter adjustments:

- Limit the RT range.
- Shorten the signal range.
- Specify a lower value for the Number of Top Matches.
- Specify a Simple Search instead of an Exhaustive Search.
- Specify Top Peaks instead of All Peaks.
- Limit the number of search types.

Figure 87. Submit Options dialog box

Submit Options

User name:

Samples:

Workflow steps to perform

☒ Acquire data 1 unacquired sample(s)

☒ Process data

Target Screening : ☒ Peak Detect ☐ Identify and Confirm

☐ Create reports

Submit as a priority sequence?

☐ Priority Sequence

☐ Next Available Batch

☒ Next Available Sample

Acquisition

Device Name	Use	Start Device
MS	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>

☒ Start when ready

Post-run system state: On

Hide Details ^ OK Cancel

Table 49. Submit Options dialog box parameters (Sheet 1 of 2)

Parameter	Description
User Name	Name of the current user.
Samples	Number of samples to be submitted for acquisition, processing, or reporting.
Workflow Steps to Perform	
Acquire Data	<p>Submits the current batch to acquisition.</p> <ul style="list-style-type: none"> When all submitted samples have been previously acquired, this option is (by default) not selected. When one or more samples in the batch have not been acquired, this option is (by default) selected.
Process Data	<p>Processes the data for the current batch using any of the following options:</p> <p>Peak Detect: Performs peak detection. You can process the data with or without performing peak detection. For example, you might want to turn off peak detection when reprocessing samples.</p> <p>Identify and Confirm: Performs both identification and confirmation.</p> <p>Note Target Screening methods can include unknown screening features. For specifying processing data for batches that also use unknown screening features, see To submit samples in the batch.</p>
Create Reports	Creates reports for the current batch.
Submit as a Priority Sequence?	
Priority Sequence	<p>With multiplexing activated, places the batch immediately after the currently acquiring batch.</p> <p>Without multiplexing activated, specifies one of the following priority options to place the batch in the queue:</p> <p>Next Available Batch: Places the batch immediately after the currently acquiring batch.</p> <p>Next Available Sample: Places the batch immediately after the currently acquiring sample.</p> <p>Note When you select Full Sequence Submission in the Configuration console, these options are unavailable because the current batch and the current sample are, in effect, the same thing.</p>

Table 49. Submit Options dialog box parameters (Sheet 2 of 2)

Parameter	Description
Acquisition pane	
Device Name	Lists all configured instruments. If the instrument that you want to use is not configured, close the TraceFinder application, configure the instrument, and then reopen the application. You cannot configure an instrument while the TraceFinder application is running. Available only when you select the Acquire Data check box.
Use	Specifies the instruments used for this acquisition. Available only when you select the Acquire Data check box.
Start Device	Specifies the instrument that initiates the communication with the other instruments. This is usually the autosampler. Available only when you select the Acquire Data check box.
Start When Ready	Starts the specified device when all the instruments are ready to acquire data. When this is cleared, individual instruments can start at different times and then must wait for the last instrument to be ready.
Post-run System State	Specifies the system state after it acquires the last batch: On (default), Standby, or Off.
Buttons	
Hide/Show Details	Collapses or expands the acquisition details of the Submit Options dialog box.
OK	Begins the selected processes.
Cancel	Closes the Submit Options dialog box without submitting any tasks.

❖ **To view the output files**

Locate the files to view from the following directories:

The application writes saved batches to the project folder:

...\\TraceFinderData\\4.0\\Projects\\...

For each acquired sample, the application writes an RSX file to the batch Data folder:

...\\TraceFinderData\\4.0\\Projects\\...\\Data

The application saves method information to the batch Methods folder:

...\\TraceFinderData\\4.0\\Projects\\...\\Methods

The application writes the reports to the batch Reports folder:

...\\TraceFinderData\\4.0\\Projects\\...*batch*\\Reports

Saving a Batch to a New Location

You can move the current batch to a different project folder, or you can make a copy of the current batch and save the copy to a different project folder.

Follow these procedures:

- [To save a batch to another project folder](#)
- [To move a batch to another folder](#)
- [To create a new project folder](#)
- [To delete project folders](#)
- [To rename project folders](#)

❖ To save a batch to another project folder

1. Choose **File > Save > Save Batch As** from the Batch View main menu in the Analysis mode.

The [Save Batch As Dialog Box](#) opens.

2. Select a storage location.

The default storage location is C:\TraceFinderData\4.0\Projects.

3. Select or create a project folder.
4. Type a name for the new batch.

If you are saving the batch to a different folder, you must give it a unique name. You cannot overwrite an existing batch in a folder.

5. Click **Save**.

When you save the batch to a different folder, the reports reflect the original project folders and the application does not save the calibration history.

❖ To move a batch to another folder

1. Choose **File > Save > Move Batch** from the Batch View main menu in the Analysis mode.

The [Save Batch As Dialog Box](#) opens.

2. Select a storage location.

The default storage location is C:\TraceFinderData\4.0\Projects.


3. Select or create a project folder.
4. Type a name for the new batch.

You must give the batch a unique name in the new subproject folder. You cannot overwrite an existing batch.

5. Click **Save**.

When you move the batch, the reports reflect the original project and subproject folders and the application does not save the calibration history.

❖ To create a new project folder

1. In the Save Batch As dialog box, select the folder for which you will create a new lower-level folder.
 - You can select the main TraceFinderData\4.0\Projects folder and create a new folder under it.
 - You can select one of the existing folders and create a lower-level folder under it.
2. Click the **Create Folder** icon, .


The application adds a new lower-level folder to the selected folder.

3. Select the new folder name and type a name for the folder.

Folder names can contain spaces and special characters, except for the following special characters: \ / : + ? " < >

Note After you add a lower-level folder, you cannot rename the parent folder.


❖ To delete project folders

1. In the Save Batch As dialog box, select the folder to delete.
2. Click the **Delete Folder** icon, .

With no confirmation prompt, the application immediately removes the selected folder.

Note This feature is not available for folders that contain lower-level project or batch folders; you must first delete the lower-level project or batch folders.

❖ To rename project folders

1. In the Save Batch As dialog box, select the folder to rename.
2. Click the **Rename Folder** icon, .

Note This feature is not available for folders that contain lower-level project or batch folders; you must first delete the lower-level project or batch folders.

3. Type a new name for the folder and press ENTER.

The application saves the new folder name.

5 Using the Analysis Mode for Target Screening Batches

Working in the Batch View for Target Screening Batches

Save Batch As Dialog Box

Use the features in the Save Batch As dialog box to save a batch to a new name or to move a batch to a different project folder.

Figure 88. Save Batch As dialog box

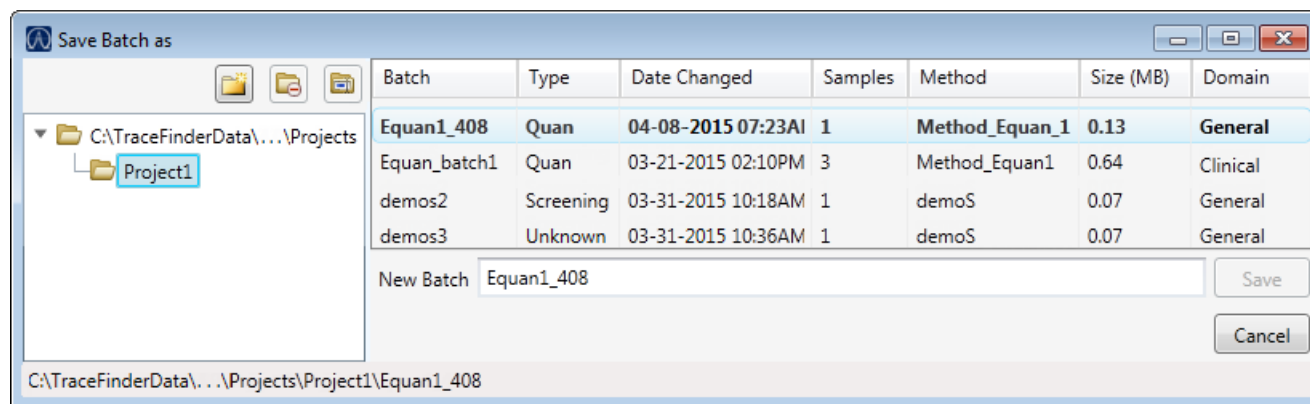


Table 50. Save Batch As dialog box parameters (Sheet 1 of 2)




Parameter	Description
 Create New Folder	Adds one of the following: <ul style="list-style-type: none">When a drive is selected, adds a new project-level folder to the drive.When a project folder is selected, adds a subproject-level folder to the selected project.When a subproject folder is selected, adds a lower-level folder to the subproject. Or, you can right-click and choose Create Folder.
 Delete Folder	With no confirmation prompt, immediately removes the selected folder. You cannot delete a folder that contains lower-level project or batch folders; you must first delete the lower-level project or batch folders. Or, you can right-click and choose Delete.
 Rename Folder	Renames the selected folder. You cannot rename a folder that contains lower-level project or batch folders; you must first delete the lower-level project or batch folders. Or, you can right-click and choose Rename.
Batch table	
Batch	Name of batches in the selected project.
Type	Type of batch: Quan, Screening, or Unknown Only.
Date Changed	Date that the batch was last updated.
Samples	Number of samples in the batch.
Method	Name of the method used to create the batch.

Table 50. Save Batch As dialog box parameters (Sheet 2 of 2)

Parameter	Description
Size	Size of the batch in megabytes.
Domain	TraceFinder domain in which the batch was created.
New batch parameters	
New Batch	Name of the new batch to create. Note If the Create button is not activated, you have entered a name that is already used or you have not selected a method.
<i>Path</i>	Path to the project in the TraceFinderData\4.0\Projects folder where the batch is created.
Buttons	
Save	Saves the batch to the specified name and folder and opens the Batch View for the new batch.
Cancel	Closes the Save Batch As dialog box without saving the batch.
Shortcut menu commands	
Create Folder	Adds one of the following: <ul style="list-style-type: none"> • When a drive is selected, adds a new project-level folder to the drive. • When a project folder is selected, adds a subproject-level folder to the selected project. • When a subproject folder is selected, adds a lower-level folder to the subproject.
Delete Folder	Immediately removes the selected folder. There is no prompt to confirm that you want to delete the selected folder. You cannot delete a folder that contains lower-level project or batch folders; you must first delete the lower-level project or batch folders.
Rename Folder	Renames the selected folder. You cannot rename a folder that contains lower-level project or batch folders; you must first delete the lower-level project or batch folders.
Expand Child Nodes	Expands all project and subproject folders in the Project tree.
Collapse Child Nodes	Collapses all project and subproject folders in the Project tree.

Working in Data Review for Target Screening Batches

The Data Review view displays the data generated by the master method. Use Data Review to verify the data for a compound before you generate reports.

Follow these procedures:

- [To open the Data Review page](#)
- [To display or hide a pane](#)
- [To move, dock, or float Data Review panes](#)
- [To restore the default layout](#)

❖ To open the Data Review page

1. Click **Analysis** in the navigation pane.

The Analysis navigation pane opens.

2. Click **Data Review**.

The Data Review view opens.

On the Data Review page, the panes show the following:

- A list of all samples in the current batch
- The compound results for all compounds in the method
- Chromatogram and spectrum plots for all compounds found in the currently selected sample

❖ To display or hide a pane

From the View menu, choose to display—indicated with a check mark—or hide the following panes:

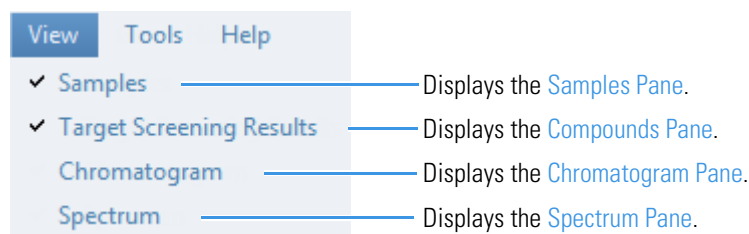
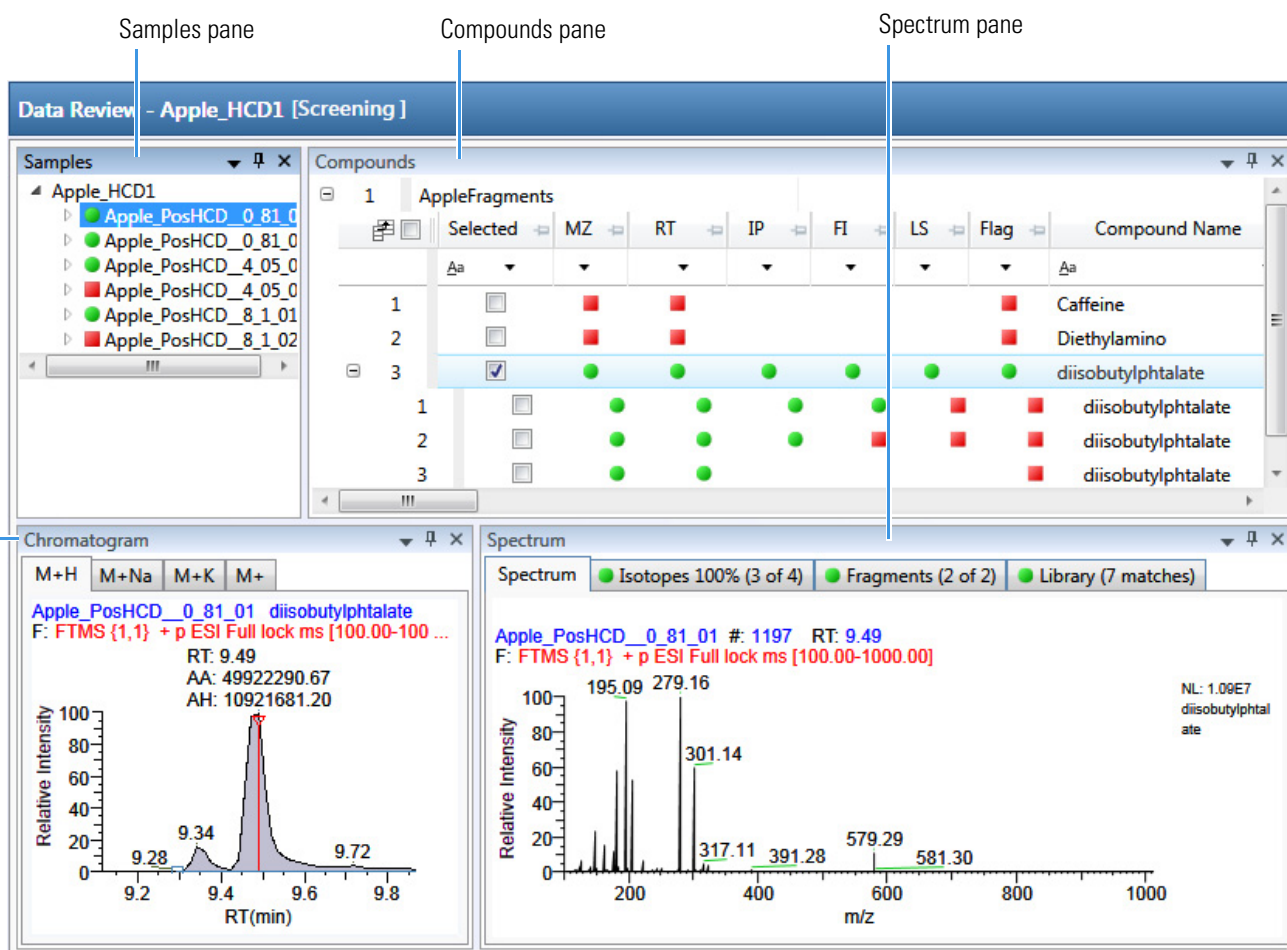


Figure 89. Data Review panes

Chromatogram pane



When you select a sample in the Samples pane, the associated Compounds pane flags any compound with errors in the selected sample. The associated Chromatogram pane displays the chromatogram, retention time, area, height, and signal-to-noise ratio for all compounds in the selected sample. The Spectrum pane highlights the compound selected in the Compounds pane. You can display, hide, or move any of these panes.

❖ **To move, dock, or float Data Review panes**

Follow the instructions in [Appendix C, “Moving Data Review Panes.”](#)

❖ **To restore the default layout**

Choose **View > Restore Default Layout**.

The Target Screening view displays the default panes in their default locations.

The Target Screening view includes the following features:

- [Samples Pane](#)
- [Compounds Pane](#)
- [Chromatogram Pane](#)
- [Spectrum Pane](#)

Samples Pane

Use the Samples pane to select a specific sample in the batch. The associated [Compounds Pane](#) displays all compounds in the method and flags any compound with errors in the selected sample.

Flags in the Samples pane indicate one of the following:

- A green circle means that the sample/compound/peak combination is identified and fully confirmed.
- ▲ A yellow triangle means that the sample/compound/peak combination is identified but not fully confirmed.
- A red square means that the sample/compound/peak combination is not identified.

Figure 90. Samples pane

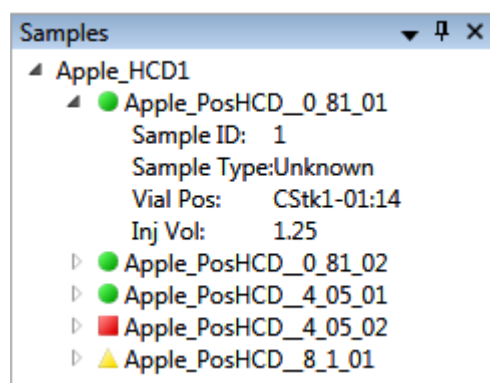


Table 51. Samples pane shortcut menu commands

Command	Description
Sort by Alphabetical	Sorts the samples alphabetically by sample name.
Sort by Import Order	Sorts the samples in the order they were processed.

Compounds Pane

The Compounds pane displays all found peaks in the selected sample and flags any compound with errors. The compounds table reflects the identified compounds found in the compound database and the results of the method processing criteria. See [Compounds Pane Columns](#).

Use the following Compounds pane features:

- [Color Coding for Measured or Calculated Values](#)
- [Displaying Multiple Adducts](#)
- [Exporting Compounds](#)
- [Compounds Pane Columns](#)

Color Coding for Measured or Calculated Values

The Compounds pane uses color-coded text to indicate the following:

- Green—Indicates that the measured value of scoring and confirmations pass the criteria specified in the method.
- Red—Indicates that the measured or calculated value does not pass the criteria specified in the method.

Measured Area	Library Score (%)
3.1312E04	12
1.9131E05	10
6.7147E04	10
1.4341E05	11

Displaying Multiple Adducts

When the application finds multiple adducts at the same retention time in a sample, the Compounds pane displays the adducts on separate rows in the table.

3	<input checked="" type="checkbox"/>	●	●	●	●	●	●	diisobutylphtalate	diisobutylphtalate@RT 9.49	C16H22O4	M+H
1	<input type="checkbox"/>	●	●	●	●	■	■	diisobutylphtalate	diisobutylphtalate@RT 9.49	C16H22O4	M+Na
2	<input type="checkbox"/>	●	●	●	■	■	■	diisobutylphtalate	diisobutylphtalate@RT 9.49	C16H22O4	M+K
3	<input type="checkbox"/>	●	●					diisobutylphtalate	diisobutylphtalate@RT 9.49	C16H22O4	M+

Exporting Compounds

Use the commands in the File menu to export data to an Excel spreadsheet, a CSV file, a compound database, or a new quantitation method.

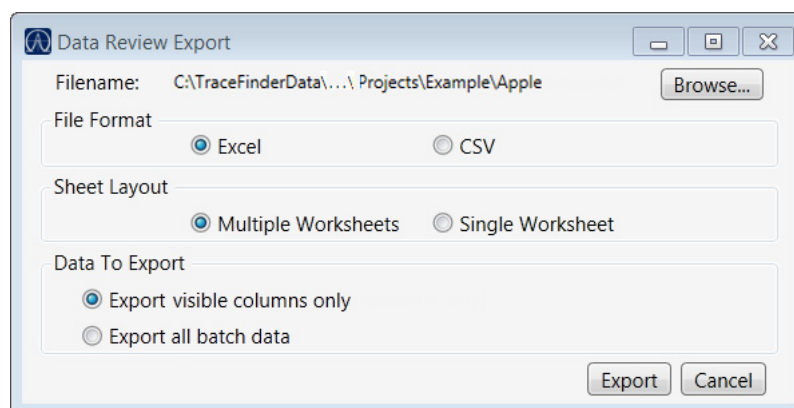
Follow these procedures:

- [To export compounds to an Excel spreadsheet](#)
- [To export compounds to a CSV file](#)
- [To export compounds to a compound database](#)
- [To create a new quantitation method with the selected compounds](#)
- [To create a new quantitation method and update the compound database](#)

❖ To export compounds to an Excel spreadsheet

1. For each compound that you want to export to an Excel spreadsheet, select the check box in the Selected column.
2. Choose **File > Export Data To > CSV or Excel** from the main menu

The Data Review Export dialog box opens.



3. Click **Browse** and, in the Export Data to Excel dialog box, locate the folder where you want to save the file.
4. Type a file name for the XLSX file and click **Save**.
5. In the File Format area, select the **Excel** option.
6. In the Sheet Layout area, select one of the following file formats for the spreadsheet.
 - **Multiple Worksheets:** Writes one sample to each Excel worksheet tab.
 - **Single Worksheet:** Writes all samples to a single Excel worksheet tab.
7. In the Data to Export area, select one of the following sets of data to export.
 - **Export Visible Columns Only:** Writes data from the displayed columns of selected samples to the specified worksheet format.

- **Export All Batch Data:** Writes data from all columns (displayed or hidden) of all samples to the specified worksheet format.

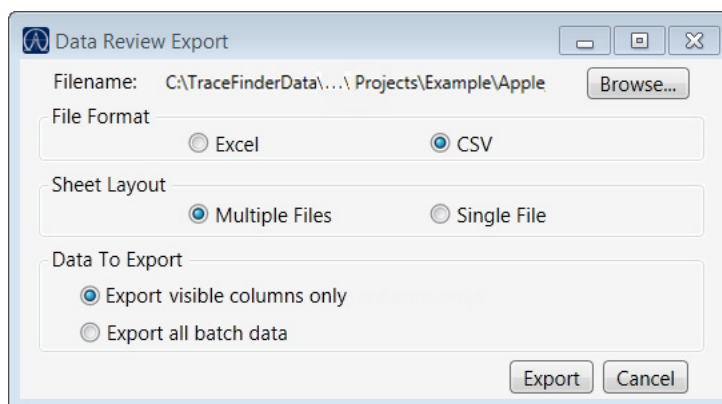
8. Click **Export**.

The application saves the specified compound data to an Excel spreadsheet and opens the folder where you saved the file.

❖ **To export compounds to a CSV file**

1. For each compound that you want to export to a CSV file, select the check box in the Selected column.
2. Choose **File > Export Data > To CSV or Excel** from the main menu.

The Data Review Export dialog box opens.



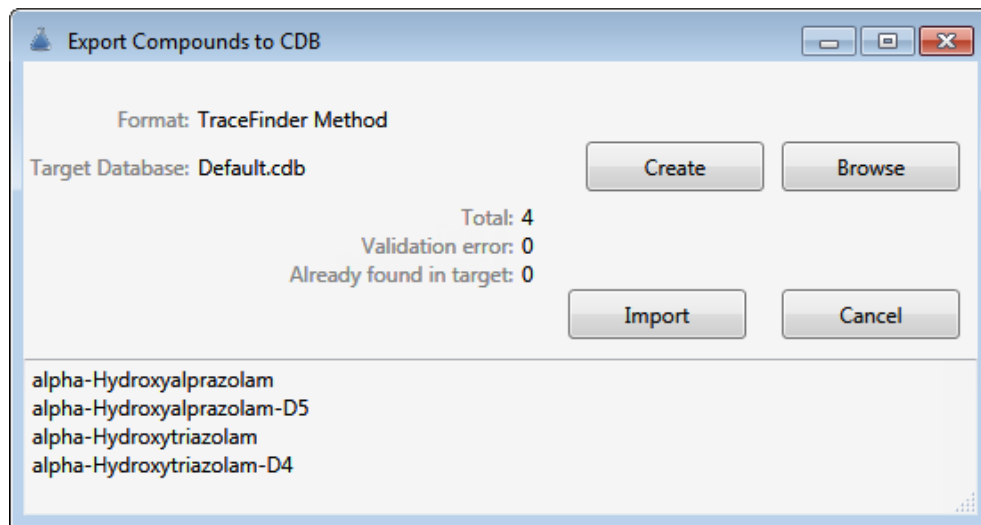
3. Click **Browse** and, in the Export Data to Excel dialog box, locate the folder where you want to save the file.
4. Type a file name for the CSV file and click **Save**.
5. In the File Format area, select the **CSV** option.
6. In the Sheet Layout area, select one of the following file formats for the spreadsheet.
 - **Multiple Files:** Writes one sample to each CSV file.
 - **Single File:** Writes all samples to a single CSV file.
7. In the Data to Export area, select one of the following sets of data to export.
 - **Export Visible Columns Only:** Writes data from the displayed columns of selected samples to the specified worksheet format.
 - **Export All Batch Data:** Writes all data from all samples in the batch to the specified worksheet format.
8. Click **Export**.

The application saves the specified compound data to an CSV spreadsheet and opens the folder where you saved the file.

❖ To export compounds to a compound database

1. For each compound that you want to export to a compound database, select the check box in the Selected column, including any adducts that you want to export.
2. Choose **File > Export Data To > Compound Database** from the main menu.

The Export Compounds to CDB dialog box opens.



3. Do one of the following:
 - Accept the default target database.
 - Click **Create** and type the name for a new compound database.
 - Click **Browse** and select from the list of compound databases.
4. Click **Import**.
 - When you export compounds to a database that already contains these compounds, the application updates the retention times in the database.
 - When you add compounds to a database that does not contain these compounds, the application adds all the compound data to the database.

When you export only one adduct for a compound, the application uses the selected adduct as the peak in the updated or new compound database. When you export multiple adducts for export, the application uses the adducts in the order of intensity.

The application does the following:

- Uses the measured retention time value for the compound in the updated or new compound database.
- Uses the expected m/z value for the compound in the updated or new compound database.
- Exports all found fragments to the updated or new compound database.

❖ **To create a new quantitation method with the selected compounds**

1. For each compound that you want to export to a new quantitation method, select the check box in the Selected column including any adducts that you want to export.
2. Choose **File > Export Data To > New Quantitation Method** from the main menu.

The Acquisition page of a new quantitation method opens.

3. From the Instrument Method list, select a method (.meth) file to use for acquiring the data.
4. Choose **File > Save** from the main menu.
5. In the Save Master Method dialog box, type a name for the method and click **OK**.
6. Click **Compounds** in the Method View navigation pane.

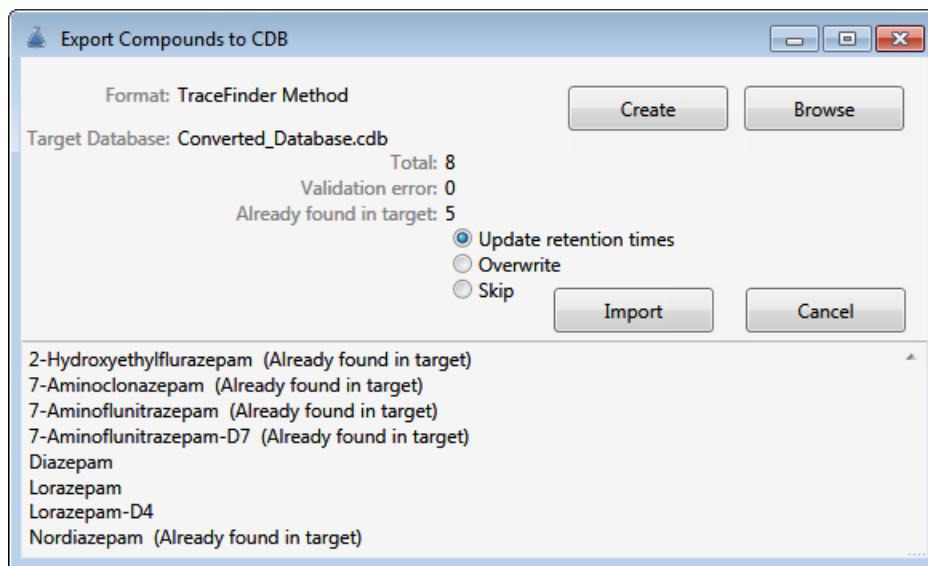
The application exports the selected compounds from the target screening method to the new quantitation method and uses data from the selected compounds as follows:

- Exports quantitation peaks in the order of intensity.
- Exports measured retention time value for the compounds in the new method.
- Exports expected m/z value for the compounds in the new method.
- Exports all found fragments to the new method.
- Adds a filter for both quantitative peaks and confirming ions.

❖ To create a new quantitation method and update the compound database

1. For each compound that you want to export, select the check box in the Selected column, including any adducts that you want to export.
2. Choose **File > Export Data > Update Compound Database and Create New Quantitation Method** from the main menu.

The Import Compounds to CDB dialog box opens.



The dialog box lists all compounds selected in the screening batch.

3. Do one of the following:
 - Accept the default target database.
 - Click **Create** and type the name for a new target database.
 - Click **Browse** and select from the list of compound databases.

The compounds list indicates which compounds already exist in the target compound database and which do not.

4. When any of the compounds already exist in the target database, choose one of the following options:
 - **Update Retention Times:** Updates only the retention times for the duplicate compounds in the target database.
 - **Overwrite:** Overwrites all compound data for the duplicate compounds in the target database.
 - **Skip:** Does not write any data from the duplicate compounds to the target database.

5. Click **Import**.

The Acquisition page of a new quantitation method opens.

Method View - Unnamed*

Calibration file last used:

Lab Name: Default Laboratory

Assay type: Assay name

Injection Volume: 1.00

Mass Precision: 2.00

Ion range calc method: Manual

Instrument Method: [Dropdown]

Edit Update


6. From the Instrument Method list, select a method (.meth) file to use for acquiring the data.
7. Choose **File > Save** from the main menu.
8. In the Save Master Method dialog box, type a name for the method and click **OK**.
9. Click **Compounds** in the Method View navigation pane.

The application exports the selected compounds (with the specified options) from the target screening method to the new quantitation method.

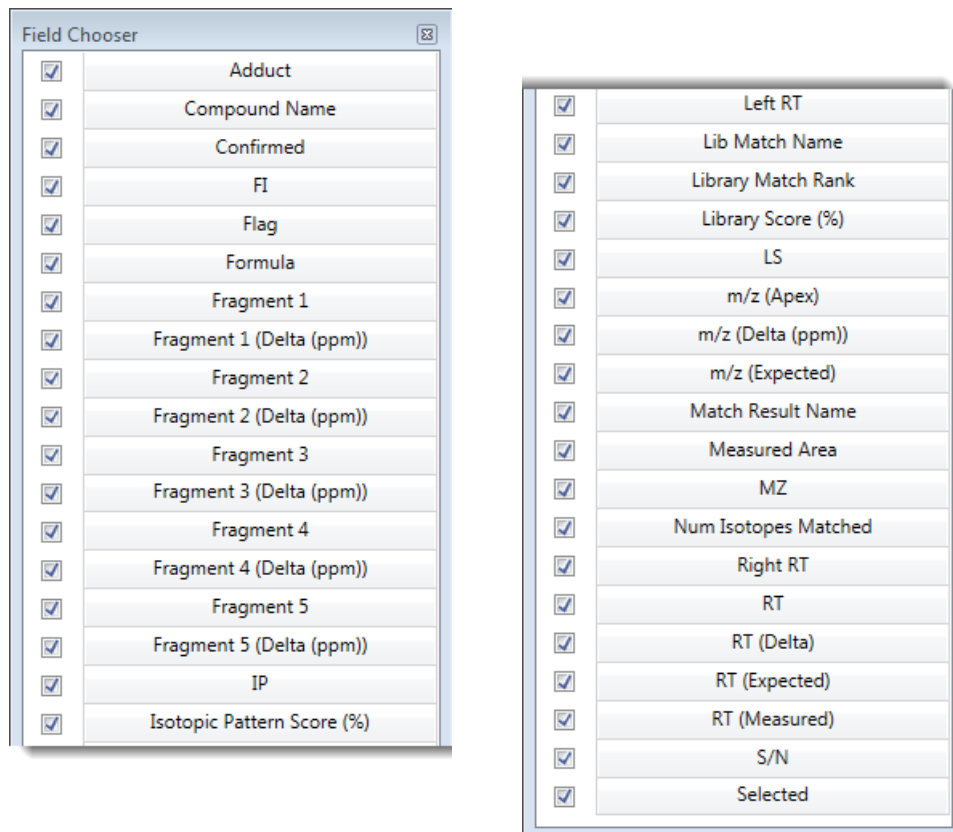
Compounds Pane Columns

The columns of data in the Compounds pane display all parameter values associated with each compound in the selected sample. See [Compounds Pane Parameters](#).

❖ To hide or display columns in the Compounds pane


1. Click the **Field Chooser** icon, , in the upper left corner of the pane.

The Field Chooser displays all available columns of data for the Compounds pane.



2. Select the check box for each column that you want to display, or clear the check box for each column that you want to hide.

The application immediately displays or hides the column in the Compounds pane.

3. When you are finished modifying the column display, click  to close the Field Chooser.

Compounds Pane Parameters

The Compounds pane displays all parameter values associated with each compound in the selected sample.

Figure 91. Compounds pane

	Selected	MZ	RT	IP	FI	LS	Flag	Compound Name	Match Result Name	Formula
1	<input type="checkbox"/>	●	●	●	●	●	●	2-Hydroxyethylflurazepam	2-Hydroxyethylflurazepam@RT 0.19	C17H14ClFN2O2
2	<input type="checkbox"/>	●	●	●	●	●	●	2-Hydroxyethylflurazepam	2-Hydroxyethylflurazepam@RT 1.42	C17H14ClFN2O2
3	<input type="checkbox"/>	●	■	●	■	■	▲	2-Hydroxyethylflurazepam	2-Hydroxyethylflurazepam@RT 1.71	C17H14ClFN2O2
4	<input type="checkbox"/>	●	■	●	■	■	▲	2-Hydroxyethylflurazepam	2-Hydroxyethylflurazepam@RT 1.96	C17H14ClFN2O2
5	<input type="checkbox"/>	●	●	●	●	●	●	2-Hydroxyethylflurazepam	2-Hydroxyethylflurazepam@RT 2.35	C17H14ClFN2O2

Table 52. Compounds pane parameters (Sheet 1 of 5)

Column	Description
	Displays the current compound database. When you have multiple screening databases, the application lists the results for each database separately.
	Expand the list of found compounds in the screening database.
Adduct	The most intense adduct for the retention time for a compound.
Compound Name	The compound name match in the compound database.
Confirmed	The number of criteria confirmed out of the total number specified in the method.
FI	<p>Fragment Ions flag. The application displays one of these indicators:</p> <ul style="list-style-type: none"> A green circle (pass) when the measured m/z value of any of the fragments is within the mass tolerance specified in the method. On the Isotopes page in the Spectrum pane, the All Isotopes and Multi-Isotopes flags are also green. A red square (fail) when the measured m/z value of none of the fragments is within the mass tolerance specified in the method. On the Isotopes page in the Spectrum pane, the All Isotopes and Multi-Isotopes flags are also red. A blank when there are no fragments detected. <p>To display a list of fragments and their pass/fail status, hold your cursor over the indicator.</p>

Table 52. Compounds pane parameters (Sheet 2 of 5)

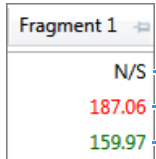
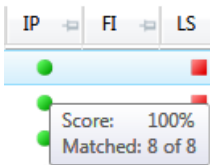
Column	Description
Flag	<p>Indicates the status of the identification and confirmation criteria.</p> <ul style="list-style-type: none"> A green circle when the sample/compound/peak combination is identified and fully confirmed. A yellow triangle when the sample/compound/peak combination is identified but not fully confirmed. A red square when the sample/compound/peak combination is not identified.
Formula	The formula for the peak as specified in the compound database.
Fragment <i>n</i>	<p>Displays the measured <i>m/z</i> for the fragment ion. The application displays a separate column for each found fragment.</p> <ul style="list-style-type: none"> For each fragment found in the compound database that passes the filter in the method, the Compounds table displays the <i>m/z</i> value in green text. For each fragment found in the compound database that does not pass the filter in the method, the Compounds table displays the <i>m/z</i> value in red text. For each fragment that is not found in the compound database, the Compounds table displays N/S (none specified). <div data-bbox="479 892 1214 1050">  <p>Fragment 1</p> <p>N/S — Fragment not found in the compound database</p> <p>187.06 — Fragment found but does not meet method parameters</p> <p>159.97 — Fragment found and meets method parameters</p> </div> <p>Note Compounds can have a maximum of five fragments, and the Compounds table has a maximum of five Fragment columns. When a compound contains fewer than five fragments, all remaining Fragment columns display N/S.</p>
Fragment <i>n</i> (Delta (ppm/mmu))	<p>The difference between the expected fragment ion <i>m/z</i> from the compound database and the measured fragment ion <i>m/z</i>.</p> <p>The application displays a separate delta column for each identified fragment.</p>
IP	<p>Isotopic Pattern flag. The application displays one of these indicators:</p> <ul style="list-style-type: none"> A green circle (pass) when the score percentage is higher than the specified fit threshold percentage. A red square (fail) when the score percentage is lower than the specified fit threshold percentage. A blank when the parameter is not scored. <p>To display the score of matched isotopes, hold your cursor over the indicator.</p> <div data-bbox="479 1617 690 1780">  <p>IP FI LS</p> <p>Score: 100%</p> <p>Matched: 8 of 8</p> </div>
Isotopic Pattern Score (%)	The percentage of the number of total isotopes to the number of matched isotopes.
Left RT	The time point of the left leading edge of the integrated peak.

Table 52. Compounds pane parameters (Sheet 3 of 5)

Column	Description
Lib Match Name	<p>The name of the best matching compound in the library search. When the application finds a match in the library, this column displays the matching library entry with the highest score.</p> <ul style="list-style-type: none"> When the application does not perform a library search, this column displays “N/A” in black text. When the application does not perform an MS/MS scan, this column displays “N/A” in red text.
Library Match Rank	<p>Displays the ranking of the library match. When the application finds a match in the library, this column displays the library entry’s relative rank, in the format “x of y”, where</p> <ul style="list-style-type: none"> x = the rank of the highest scoring library match. y = the total number of library matches from the list of matches for a particular adduct that contains the highest scoring match. <p>When the application performs both a library search and an MS/MS scan and both the library entry and the formula match the target compound:</p> <ul style="list-style-type: none"> A library score that is higher than or equal to the score threshold meets the criteria, and the values in this column appear in green. A library score that is lower than the score threshold fails to meet the criteria, and the values in this column appear in red. <p>When the application does not perform a library search, this column displays “N/A” in black text.</p> <p>When the application does not perform an MS/MS scan, this column displays “N/A” in red text.</p>
Library Score (%)	<p>The score from the library fit. When the application finds a match in the library, this column displays the highest score associated with the Lib Match Name parameter.</p> <p>When the application performs both a library search and an MS/MS scan and both the library entry and the formula match the target compound:</p> <ul style="list-style-type: none"> A library score that is higher than or equal to the score threshold meets the criteria, and the values in this column appear in green. A library score that is lower than the score threshold fails to meet the criteria, and the values in this column appear in red. <p>When the application does not perform a library search, this column displays “N/A” in black text.</p> <p>When the application does not perform an MS/MS scan, this column displays “N/A” in red text.</p> <p>Valid range: 1 through 100%</p>
LS	<p>Library Search flag. The application displays one of these flags:</p> <ul style="list-style-type: none"> A green circle when the library search is successful. A red square when the library search is not successful.

Table 52. Compounds pane parameters (Sheet 4 of 5)

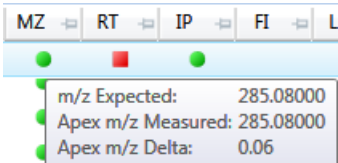
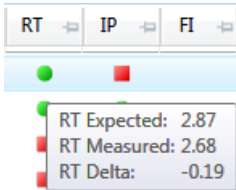
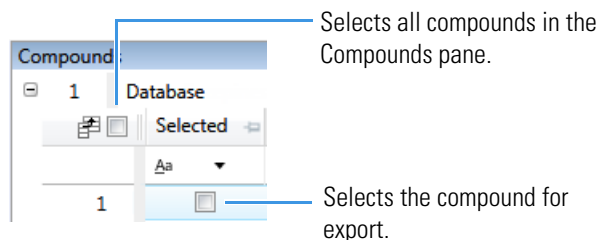
Column	Description
<i>m/z</i> (Apex)	<p>Mass-to-charge ratio found in the spectra for the peak. Assumes the charge is 1.</p> <p>When the application successfully integrates the peak, this column displays the charged <i>m/z</i> value for the compound, which is the highest intensity in the apex scan.</p> <p>When the application cannot successfully integrate the peak, this column displays “N/F” (not found).</p>
<i>m/z</i> (Delta)	<p>Difference between the <i>m/z</i> (Expected) and <i>m/z</i> (Apex). Assumes the charge is 1.</p> <p>When the <i>m/z</i> (Apex) column displays <i>m/z</i> value for the compound, this column displays the delta <i>m/z</i> corresponding to the highest intensity in the apex scan.</p> <ul style="list-style-type: none"> When the mass tolerance is specified in ppm in the master method, then $m/z \text{ (Delta)} = 1\,000\,000 \times ([m/z \text{ (Apex)} - m/z \text{ (Expected)}] \div m/z \text{ (Expected)})$. When the mass tolerance is specified in mmu in the master method, then $m/z \text{ (Delta)} = 1000 \times (m/z \text{ (Apex)} - m/z \text{ (Expected)})$.
<i>m/z</i> (Expected)	<p>Mass-to-charge ratio from the compound database. Assumes the charge is 1.</p> <ul style="list-style-type: none"> When an adduct is found, the application displays the neutral mass value for the compound (calculated from the neutral formula) \pm the mass of the most intense adduct ion found for the compound. When no adduct is found, the application displays the neutral mass value for the compound \pm the mass of the first adduct entered in the compound database. <p>For details about defining adducts for the compound database, refer to Chapter 1, “Using the Configuration Console,” in the <i>TraceFinder Lab Director User Guide</i>.</p> <p>For details about adding adducts to compounds in the database, refer to Chapter 2, “Using Compound Databases in the Method Development Mode,” in the <i>TraceFinder Lab Director User Guide</i>.</p> <p>Note When the adduct is a gain, the adduct mass is a positive number. When the adduct is a loss, the adduct mass is a negative number. The resulting mass value after adding or subtracting the adduct mass is always a positive number.</p>
Match Result Name	The compound name match in the compound database and the retention time.
Measured Area	The AA value from the chromatogram pane.
MZ	<p>Mass-to-charge ratio flag. The application displays one of these indicators:</p> <ul style="list-style-type: none"> A green circle (pass) when the measured <i>m/z</i> value is within the specified threshold. A red square (fail) when the measured <i>m/z</i> value is not within the specified threshold. A blank when the mass-to-charge value is unavailable. <p>To display the expected, measured, and delta <i>m/z</i>, hold your cursor over the indicator.</p> 

Table 52. Compounds pane parameters (Sheet 5 of 5)

Column	Description
Num Isotopes Matched	The number of isotopes matched in the expected calculated isotope spectra relative to the total number of isotopes used in the score calculation, in the format “x of y”, where <ul style="list-style-type: none"> x = the number of isotopes matching the elemental composition used for the Isotopic Pattern Score calculation. y = the total number of isotopes considered in the Isotopic Pattern Score calculation. This is the number of isotope peaks expected to be above the spectral noise.
Right RT	The time point of the right trailing edge of the integrated peak.
RT	Retention Time flag. The application displays one of these indicators: <ul style="list-style-type: none"> A green circle (pass) when the measured retention time value is within the RT Window value specified in the compound database. A red square (fail) when the measured retention time value is not within the RT Window value specified in the compound database. In turn, this results in a failure flag for the <i>m/z</i> value because the application cannot identify an <i>m/z</i> value that meets the retention time. A blank when the retention time value is unavailable. When the retention time is selected as “confirm” and the <i>m/z</i> is not detected, there is no flag. <p>Or, you can set a different retention time window on the Processing page in the method. Refer to Chapter 5, “Using the Method Development Mode for Target Screening Methods,” in the <i>TraceFinder Lab Director User Guide</i>.</p> <p>To display the expected, measured, and delta retention times, hold your cursor over the indicator.</p> 
RT (Delta)	Difference between the expected and measured retention time for the peak.
RT (Expected)	The retention time for the peak as specified in the compound database.
RT (Measured)	The found retention time for the peak apex.
S/N	The signal-to-noise ratio calculated for the found peak.
Selected	Identifies individual compounds for export. To select all compounds for export, select the check box in the first column.



Chromatogram Pane

Use the Chromatogram pane to display all extracted chromatograms of all adducts of the selected compound.

The first tab displays the most intense target adduct for the peak result. Additional (optional) tabs display extracted ion chromatograms for other adducts for the target compound at the same retention time in order of intensity. If no signal exists for an adduct, it displays the XIC of the expected m/z within the specified retention and chromatogram windows. When you do not specify a retention time or window, the application displays the full time range.

For each adduct, the [Spectrum Pane](#) displays the spectrum, isotopes, fragments, and library matches.

Figure 92. Chromatogram pane

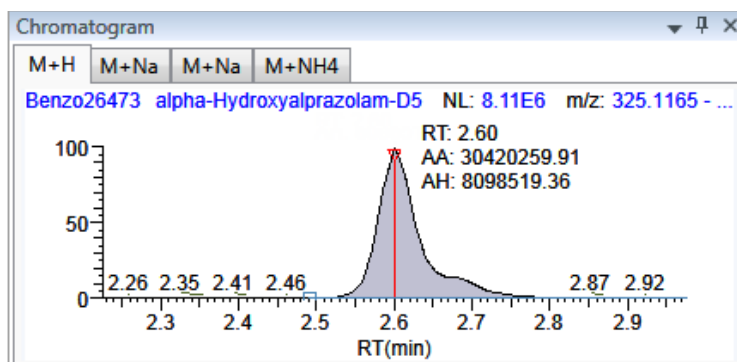


Table 53. Chromatogram pane shortcut menu commands

Command	Description
Reset Scaling	Resets the original scaling after a zoom operation.
Copy to Clipboard	Copies the graphic display to the Clipboard.

Spectrum Pane

Use the Spectrum pane to display the spectrum, isotopes, fragments, and library search information for the selected adduct in the Chromatogram pane. The Spectrum pane displays only the identification and confirmation criteria specified in the method. The confirmations are based only on the most intense adduct. Refer to “Editing the Processing Page” in Chapter 5, “Using the Method Development Mode for Target Screening Methods,” in the *TraceFinder Lab Director User Guide*.

The Spectrum pane includes the following pages of information (when available) for each selected sample/compound/peak combination:

- [Spectrum](#)
- [Isotopes](#)
- [Fragments](#)
- [Library](#)

Spectrum

The application displays the neutral loss (NL) and compound/peak name information on the right side of the [Spectrum page](#). When data is available, the plot width is the full mass range in the raw data file. Otherwise, the application scales the width to the scan range.

Follow these procedures:

- [To zoom in on a peak](#)
- [To copy a plot to another format](#)
- [To customize the labels in a small molecule or peptide spectrum](#)

❖ To zoom in on a peak

1. On the Spectrum page, drag the cursor to delineate a rectangle around the peak.
The delineated area expands to fill the view.
2. To restore the default view, right-click the Spectrum page and choose **Reset Scaling**.

❖ To copy a plot to another format

1. In the data spectrum or the reference spectrum, right-click and choose **Copy to Clipboard**.

The application copies the specified plot graphic to the Clipboard.

Note The application copies only the indicated data spectrum or the reference spectrum. It does not copy the entire Spectrum page.

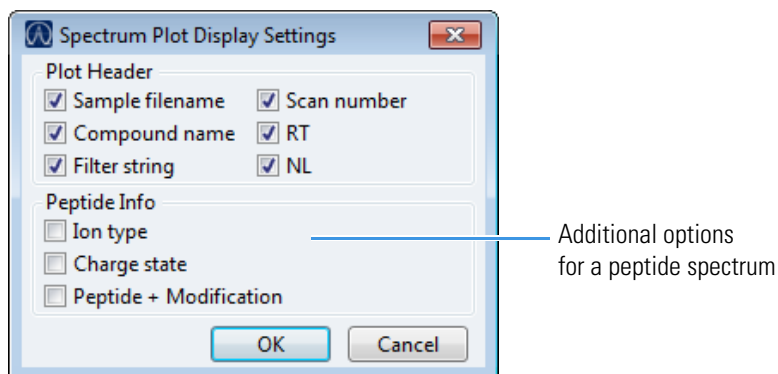
2. In the source format (for example, an Excel spreadsheet or email window), right-click and choose **Paste**.

The application pastes the graphic into the new format.

❖ To customize the labels in a small molecule or peptide spectrum

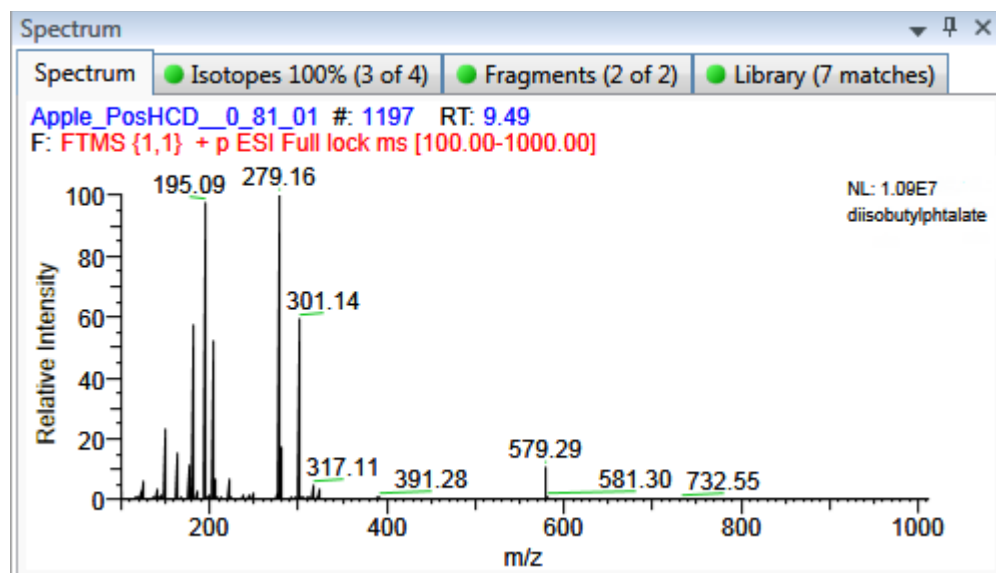
1. Right-click the [Spectrum page](#) and choose **Display Settings**.

The Spectrum Plot Display Settings dialog box opens.



2. Select the check box for each label that you want to display and click **OK**.

Figure 93. Spectrum page



Isotopes

The Isotopes page displays isotopic pattern results for all adducts of a compound according to the threshold and deviation parameters defined in the screening method.

To identify or confirm the presence of a compound, the resulting score percentage from isotopic pattern matching must be higher than the specified fit threshold percentage.

- An isotope peak is not found if its intensity, relative to the monoisotopic ion's intensity, is more than the specified intensity deviation percentage away from the theoretical relative intensity of the isotope ion.
- An isotope peak is found if its measured m/z is less than the specified mass deviation amount away from its expected m/z .

To specify threshold and deviation parameters, refer to “Editing the Processing Page” in Chapter 5, “Using the Method Development Mode for Target Screening Methods,” in the *TraceFinder Lab Director User Guide*.

The Isotopes page displays the isotopes in one of three ways:

- [All Isotopes](#)
- [Multi-Isotopes](#)
- [Individual Isotopes](#)

All isotopes pages use a shortcut menu so that you can specify how to display the data. See [Isotopes Page Shortcut Menu](#).

All Isotopes

The All Isotopes view displays a composite of all isotopes found in the compound. The application scales the window with respect to the most intense isotope. The most intense isotope is usually the first isotope unless you are using halogenated compounds. The application displays the measured peak as a solid red line; the application displays the expected peak as a dashed blue line.

The application displays these headers for the All Isotopes view:

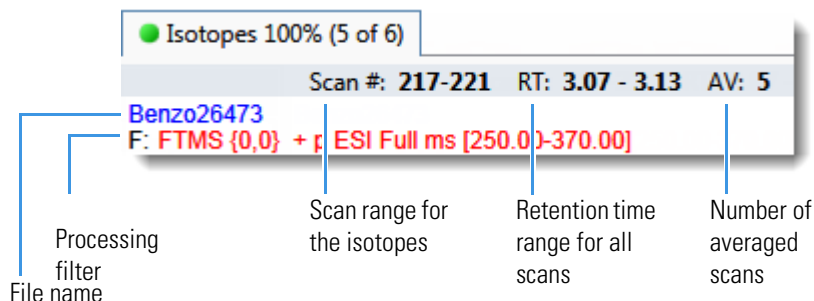


Figure 94. Isotopes page with stacked spectra for all isotopes

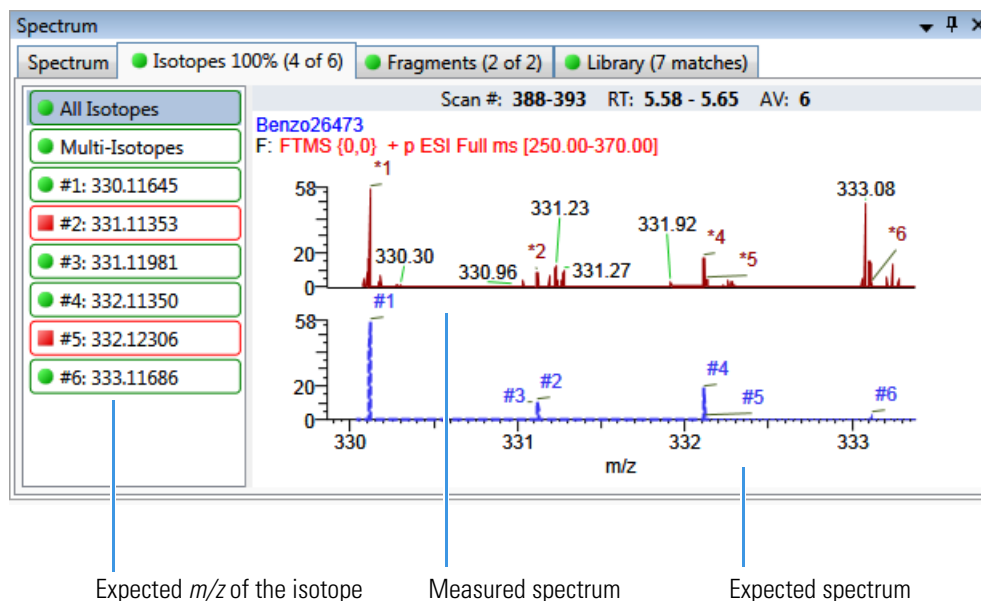
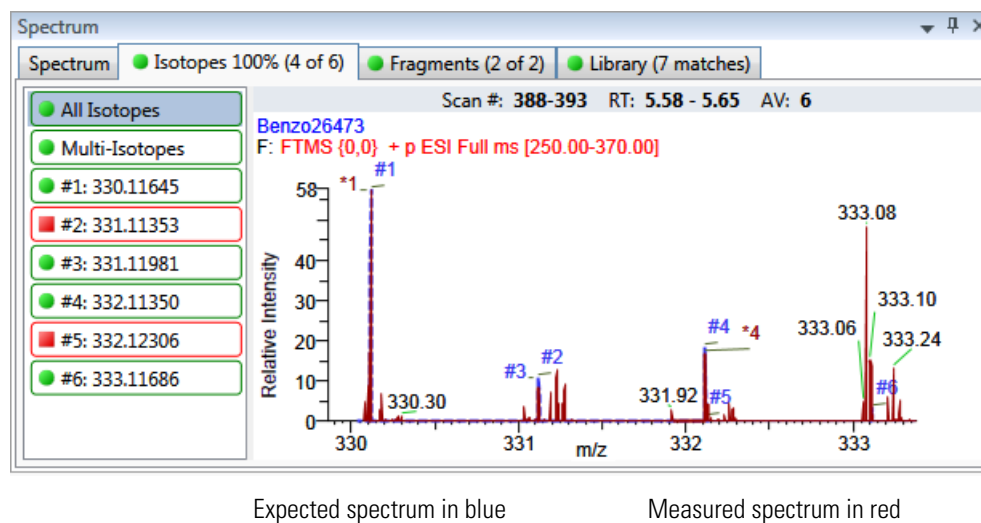


Figure 95. Isotopes page with overlaid spectra for all isotopes



Multi-Isotopes

The Multi-Isotopes view displays individual plots for each isotope. You can individually stack or overlay the plots for each isotope.

The application displays these headers for the Multi-Isotopes view:

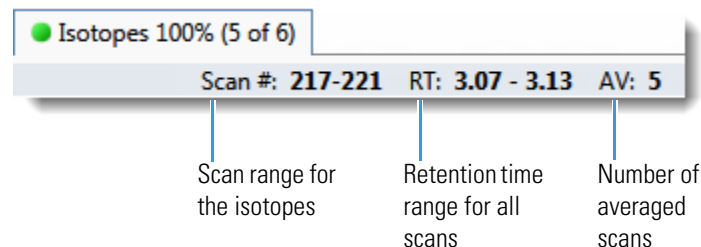


Figure 96. Isotopes page with overlaid spectra for multi-isotopes

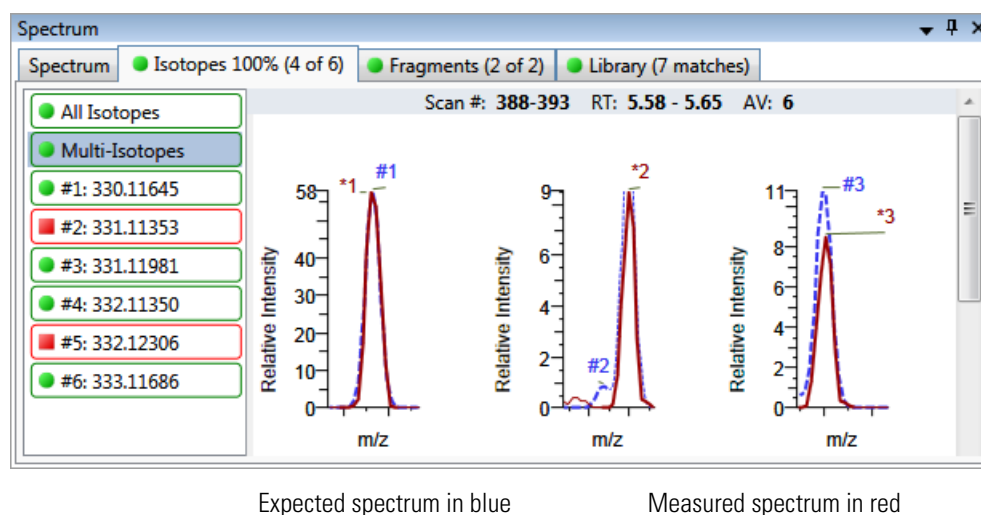
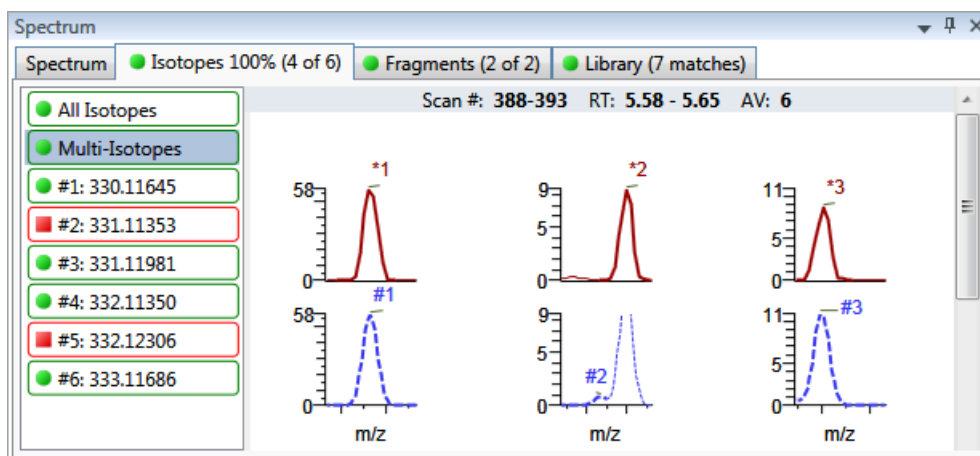


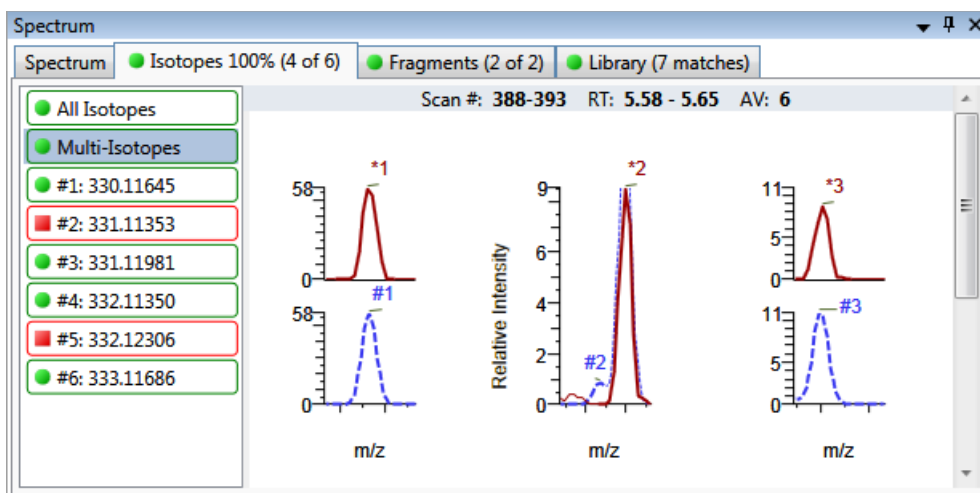
Figure 97. Isotopes page with stacked spectra for multi-isotopes



Expected spectrum in blue

Measured spectrum in red

Figure 98. Isotopes page with stacked and overlaid spectra for multi-isotopes



Expected spectrum in blue

Measured spectrum in red

Individual Isotopes

The individual isotopes view displays the expected and measured peaks for a single isotope.

The application displays these headers for the individual isotopes view:

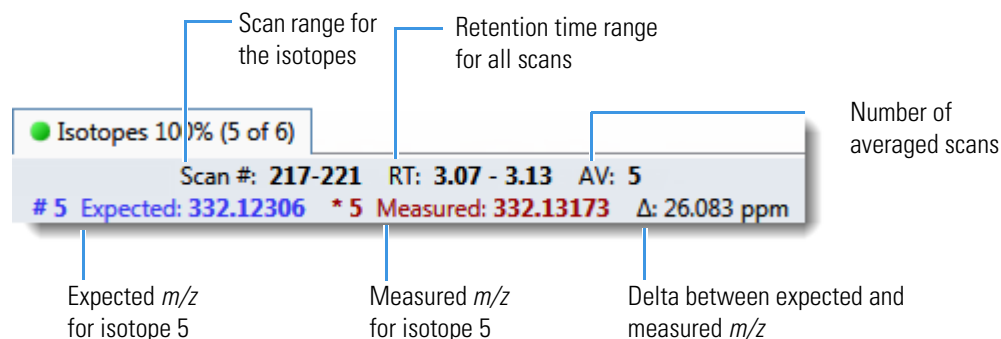


Figure 99. Isotopes page with overlaid spectra for a single isotope

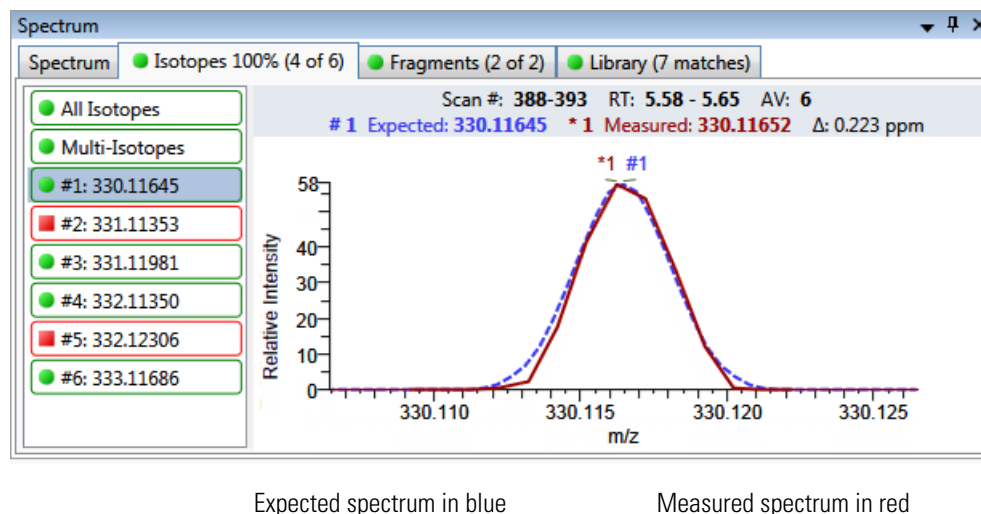
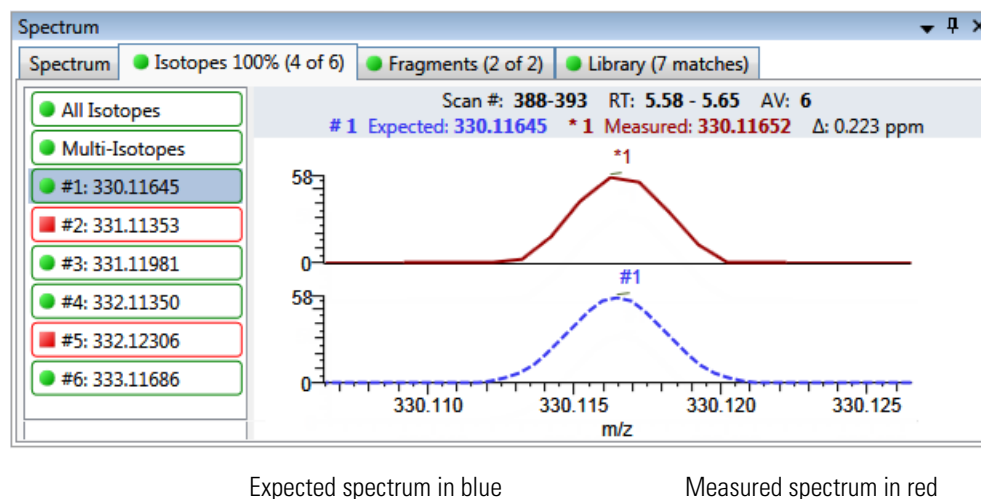


Figure 100. Isotopes page with stacked spectra for a single isotope



Isotopes Page Shortcut Menu

Use the commands in the shortcut menu to specify how you want the data displayed.

Table 54. Isotopes page shortcut menu commands

Command	Description
Reset Scaling	Resets the original scaling after a zoom operation.
Copy to Clipboard	Copies the graphic display to the Clipboard.
Display Overlay Spectra Display Stack Spectra	Overlays the two spectrum displays, or stacks the peak apex spectrum over the simulated spectrum.
Show/Hide Noise Label	Adds a noise label to each peak. Expected isotope peaks (displayed in blue) do not display a noise label.
Show/Hide Resolution Label	Adds a resolution label to each peak. Expected isotope peaks (displayed in blue) do not display a resolution label.

Fragments

The Fragments page displays the maximum number of fragments as specified in the screening method. Refer to “Editing the Processing Page” in Chapter 5, “Using the Method Development Mode for Target Screening Methods,” in the *TraceFinder Lab Director User Guide*.

If there are no fragments defined in the screening library for the compound, you can add fragments to the screening library. Refer to the instructions “To add a fragment to a target peak” in Chapter 2, “Using Compound Databases in the Method Development Mode,” in the *TraceFinder Lab Director User Guide*.

The Fragments page displays the fragments in one of two ways:

- [All Fragments](#)
- [Individual Fragments](#)

All fragments pages use a shortcut menu so that you can specify how to display the data. See [Fragments Page Shortcut Menu](#).

All Fragments

The All Fragments view displays a composite of all fragments found in the compound. The application displays the measured peak as a solid red line; the application displays the expected peak as a dashed blue line.

The application displays these headers for the All Fragments view:

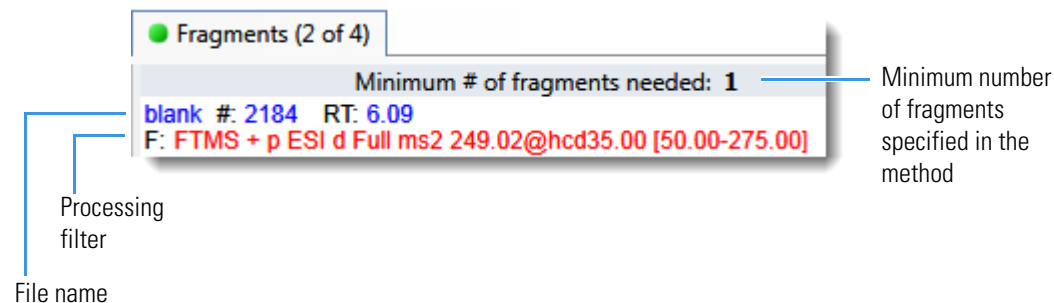


Figure 101. Fragments page with overlaid spectra for all fragments

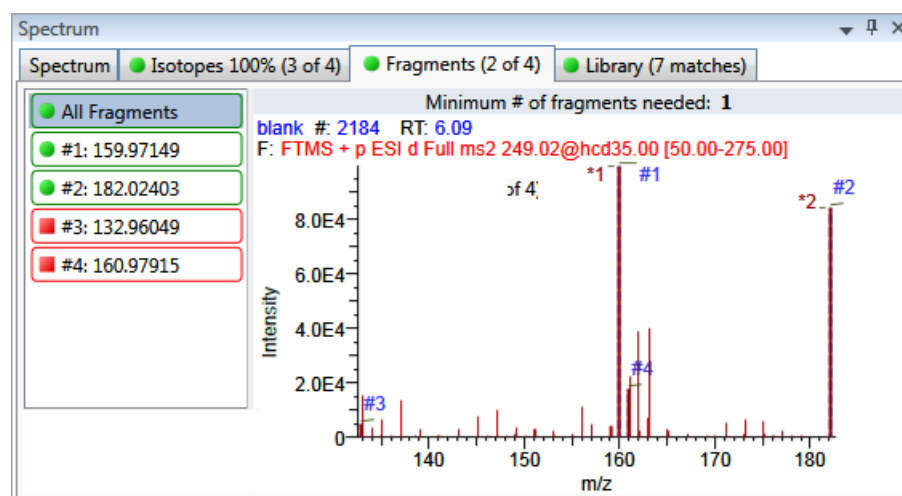
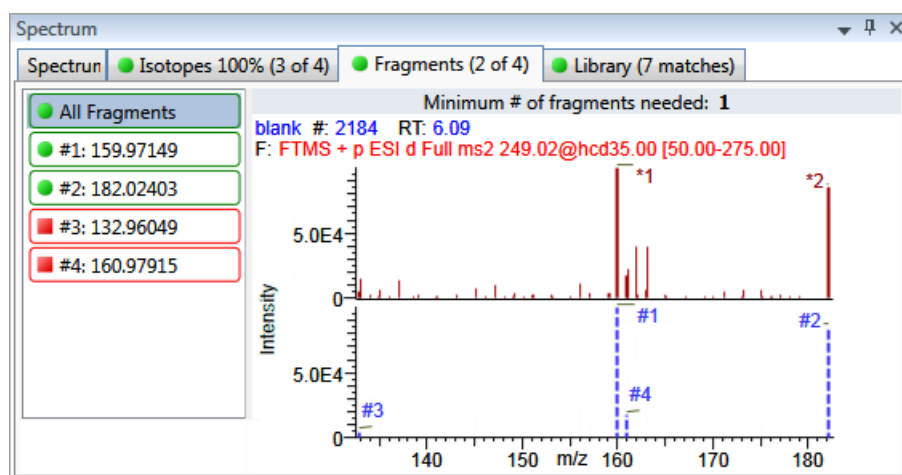


Figure 102. Fragments page with stacked spectra for all fragments



Individual Fragments

The individual fragments view displays the expected and measured peaks for a single fragment.

The application displays these headers for the individual fragments view:

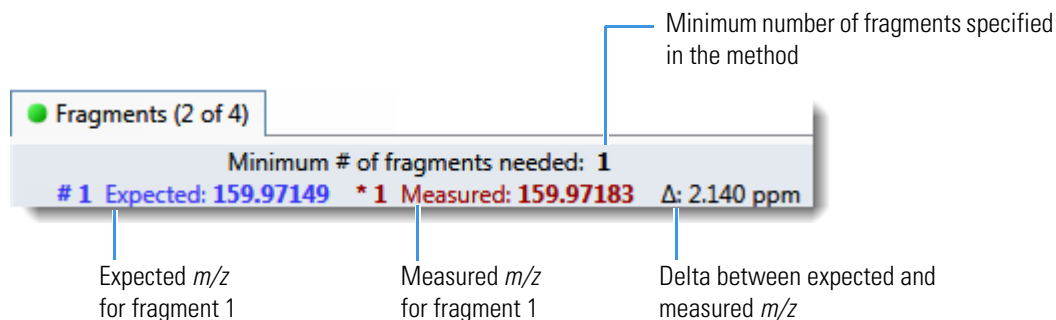


Figure 103. Fragments page with overlaid spectra for a single fragment

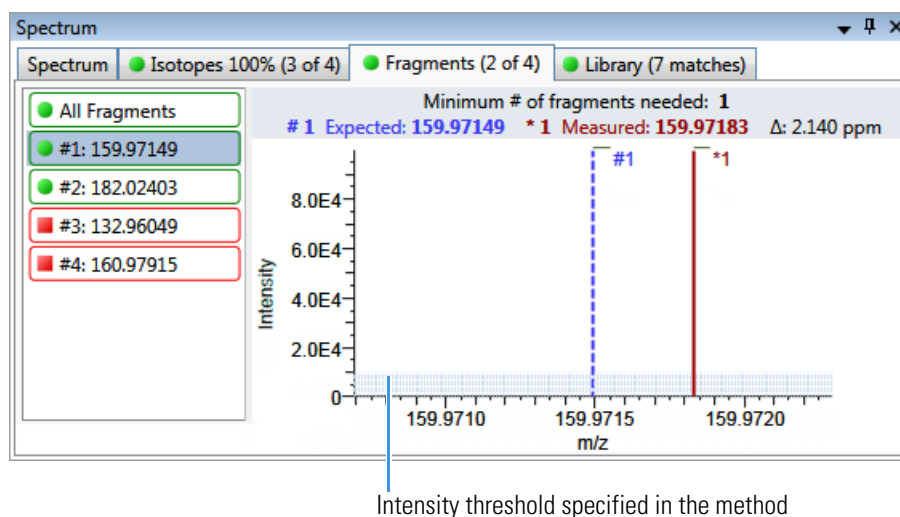
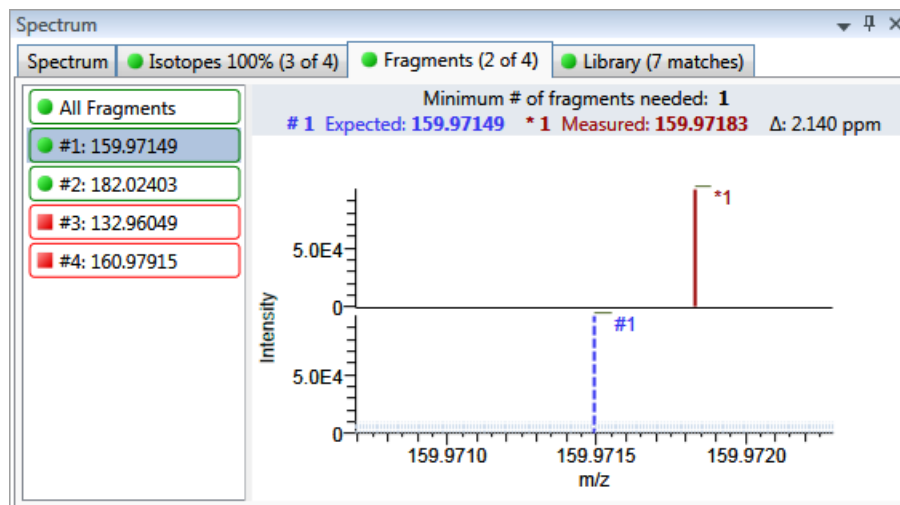


Figure 104. Fragments page with stacked spectra for a single fragment



Fragments Page Shortcut Menu

Use the commands in the shortcut menu to specify how you want the data displayed.

Table 55. Fragments page shortcut menu commands

Command	Description
Reset Scaling	Resets the original scaling after a zoom operation.
Copy to Clipboard	Copies the graphic display to the Clipboard.
Display Overlay Spectra	Overlays the two spectrum displays, or stacks the simulated spectrum and the peak apex spectrum.
Display Stack Spectra	

Library

The Library page displays the matching library spectrum (in blue) and the experimental spectrum (in black). The resulting score percentage from a library search match must be higher than your specified threshold value to identify or confirm the presence of a compound. Refer to “Editing the Processing Page” in Chapter 5, “Using the Method Development Mode for Target Screening Methods,” in the *TraceFinder Lab Director User Guide*.

The application scales both the matched library spectrum and the highest peak in the experimental spectra at 100 percent intensity and displays the resulting neutral loss (NL) value for the matched library entry name on the right side of the plot.

The application displays these headers for the individual adducts:

Library match name		Library score percentage	
● Library (7 matches)			
M+H #1: Phthalic_acid_bis-iso-butyl_ester	C16H22O4	Score: 88	Rank: 1 of 7
Adduct	Formula	Library match rank	

Figure 105. Library page with stacked spectra

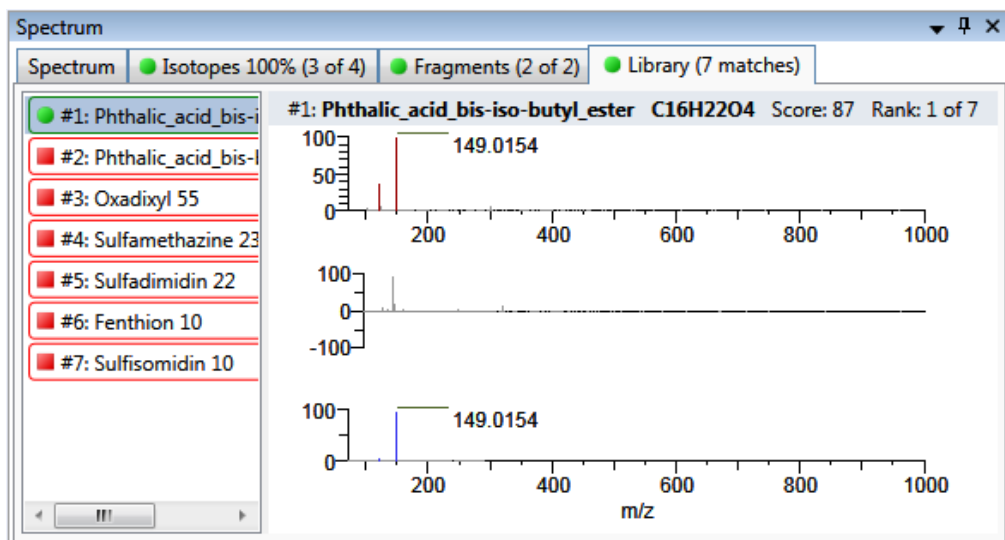
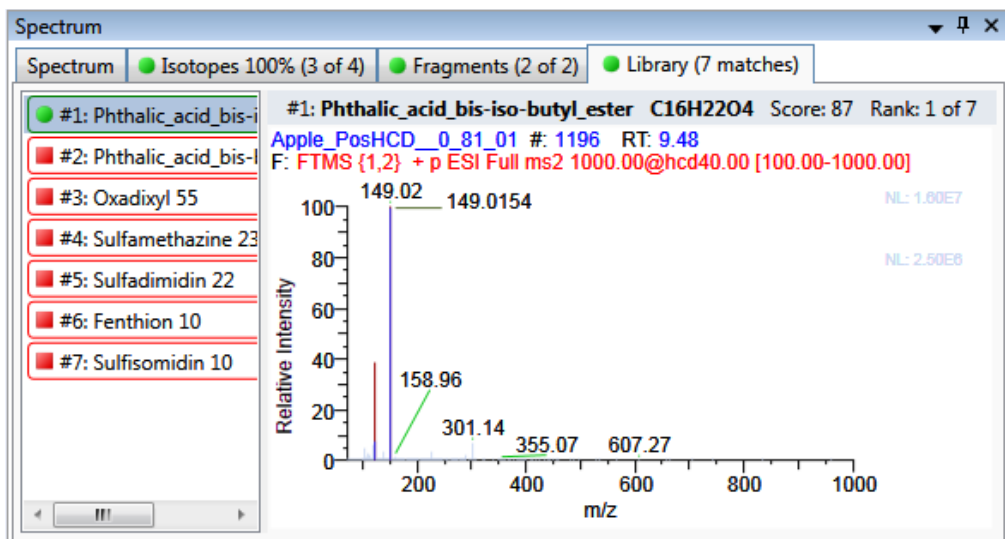


Figure 106. Library page with overlaid spectra



Library Page Shortcut Menu

Use the commands in the shortcut menu to specify how you want the data displayed.

Table 56. Library page shortcut menu commands

Command	Description
Reset Scaling	Resets the original scaling after a zoom operation.
Copy to Clipboard	Copies the graphic display to the Clipboard.
Display Overlay Spectra	Overlays the two spectrum displays, or stacks the simulated spectrum and the peak apex spectrum.
Display Stack Spectra	

Working in the Report View for Target Screening Batches

The Report View displays example reports for the current batch. You must have an open batch to use the features in the Report View.

Follow these procedures:

- [To open the Report View](#)
- [To preview a report](#)
- [To generate a report as a PDF, an Excel, or a CSV file](#)
- [To print a report](#)
- [To display a generated report](#)
- [To edit a report template](#)
- [To create a new report template](#)

❖ To open the Report View

Click **Report View** in the navigation pane.



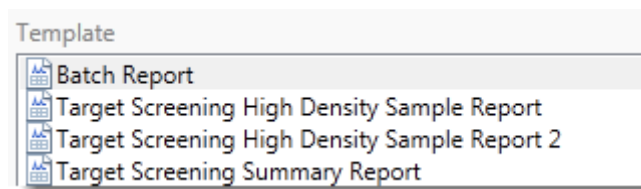
The application opens the [Report View](#).

❖ To preview a report

1. In the Template pane, select a report template.

The template list shows all the report templates that you configured in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.

Figure 107. Example of a report template list




2. Click **Preview**, .

The application opens the Report Designer, showing the report information for the current batch in the selected report template format.

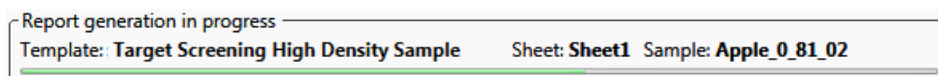
For details about using the Report Designer, see [Chapter 7, “Using the Report Designer.”](#)

❖ **To generate a report as a PDF, an Excel, or a CSV file**

1. In the Template pane, select a report template.
2. Select the check box for each of the file types that you want to create: **PDF**, **Excel**, or **CSV**.
3. Click **Generate**,  **Generate**.


The application does the following:

- Displays a green progress bar as it generates the reports.



- Creates a report for the current batch as a PDF, an Excel, or a CSV file, using the selected report template format.
- Adds information about the generated report to the Generated Reports pane.
For details about the Generated Reports pane, see [Report View](#).
- Saves the report files to the ...\\TraceFinderData\\4.0\\Projects*batch*\\ReportOutput folder.

❖ **To print a report**

1. In the Template pane, select a report template.
2. Select the check box for the **Print** file format.
3. Click **Generate**,  **Generate**.

The application does the following:

- Creates a report for the current batch using the selected report template format.
- Prints the report to your default printer.
- Adds information about the generated report to the Generated Reports pane.
For details about the Generated Reports pane, see [Report View](#).
- Saves the report files to the ...\\TraceFinderData\\4.0\\Projects*batch*\\ReportOutput folder.

❖ **To display a generated report**

In the Generated Reports pane, click **View** for the report that you want to see.

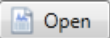
Generated Reports				
Template	Rule	Sample	Output	Generated Report File
Batch Report	Batch		pdf	View... Batch Report_20140113092939.pdf (3 pages)

Opens the generated file.

The application opens the output file.

❖ To edit a report template

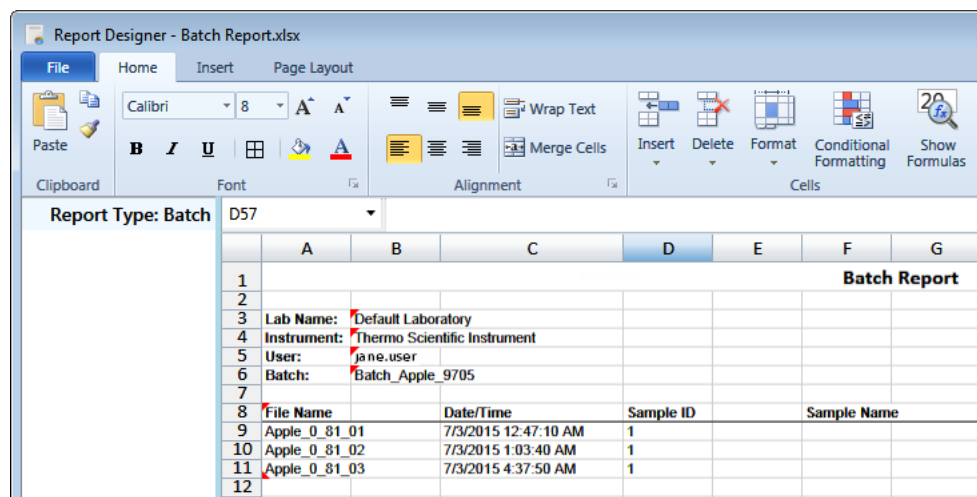
Note You cannot edit report templates that were provided with the TraceFinder application; however, you can open a TraceFinder template, make changes, and save it to a new template name.

1. In the Template pane, select a report template.
2. Click **Open**, .

The application opens the Report Designer showing the template in an Excel spreadsheet. See [Report Designer showing the template for the selected report](#).

Note When user security is activated, you must have Template Editing permission to edit report templates created by your laboratory. If the Open button is not active, user security is activated and you do not have Template Editing permission.

Figure 108. Report Designer showing the template for the selected report



3. Use the features in the Report Designer to edit the template.
For details about using the Report Designer, see [Chapter 7, “Using the Report Designer.”](#)
4. When you finish your changes, choose **File > Save** from the Report Designer menu bar.

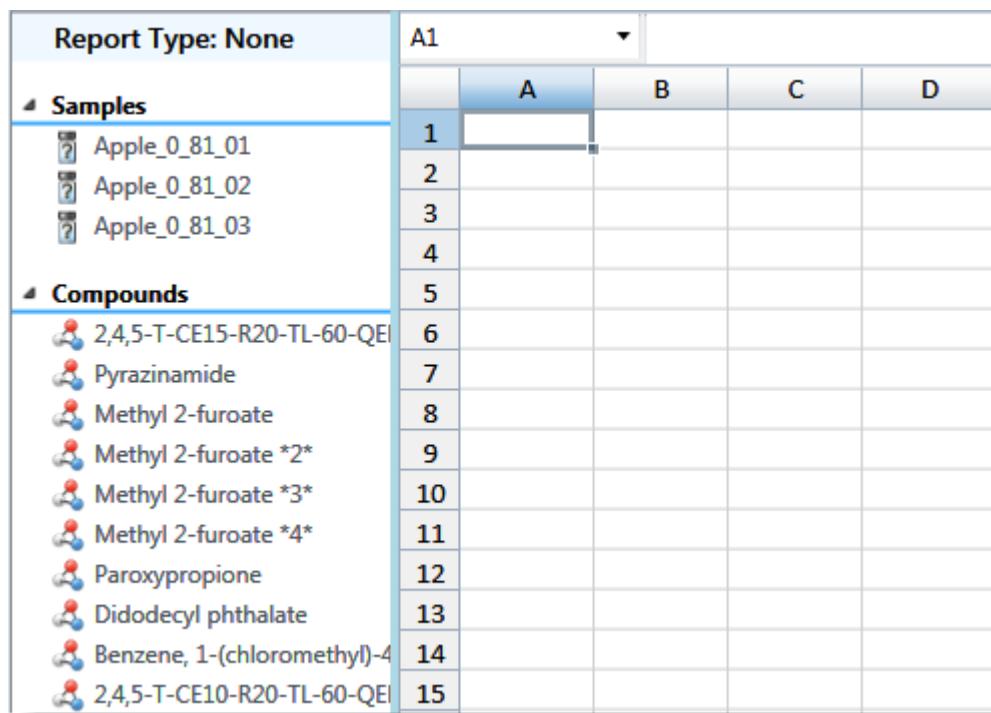
❖ To create a new report template

1. Click **New**, .

The application opens the Report Designer showing an empty template in an Excel spreadsheet. See [Report Designer showing a new, empty template](#).

The Report Type is None.

In the left pane, the spreadsheet lists all samples in the current batch and all compounds in the method used for the batch.

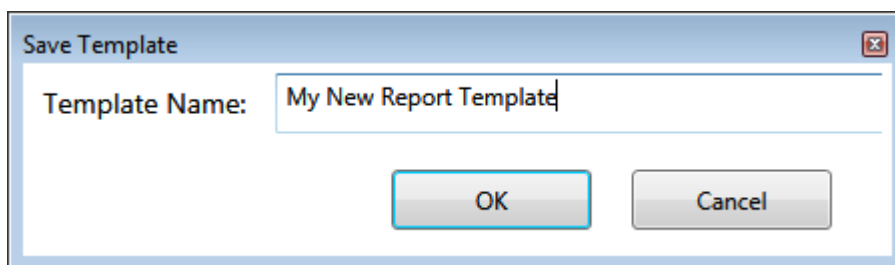
Figure 109. Report Designer showing a new, empty template

2. Use the features in the Report Designer to create the report template.

For details about using the Report Designer, see [Chapter 7, “Using the Report Designer.”](#)

3. When you finish your changes, choose **File > Save** from the Report Designer menu bar.

The Save Template dialog box opens.

Figure 110. Save Template dialog box

4. Type a name for the new report template and click **OK**.

5 Using the Analysis Mode for Target Screening Batches

Working in the Report View for Target Screening Batches

Report View

Use the features in the Report View to display example reports for the current batch.

Figure 111. Report View

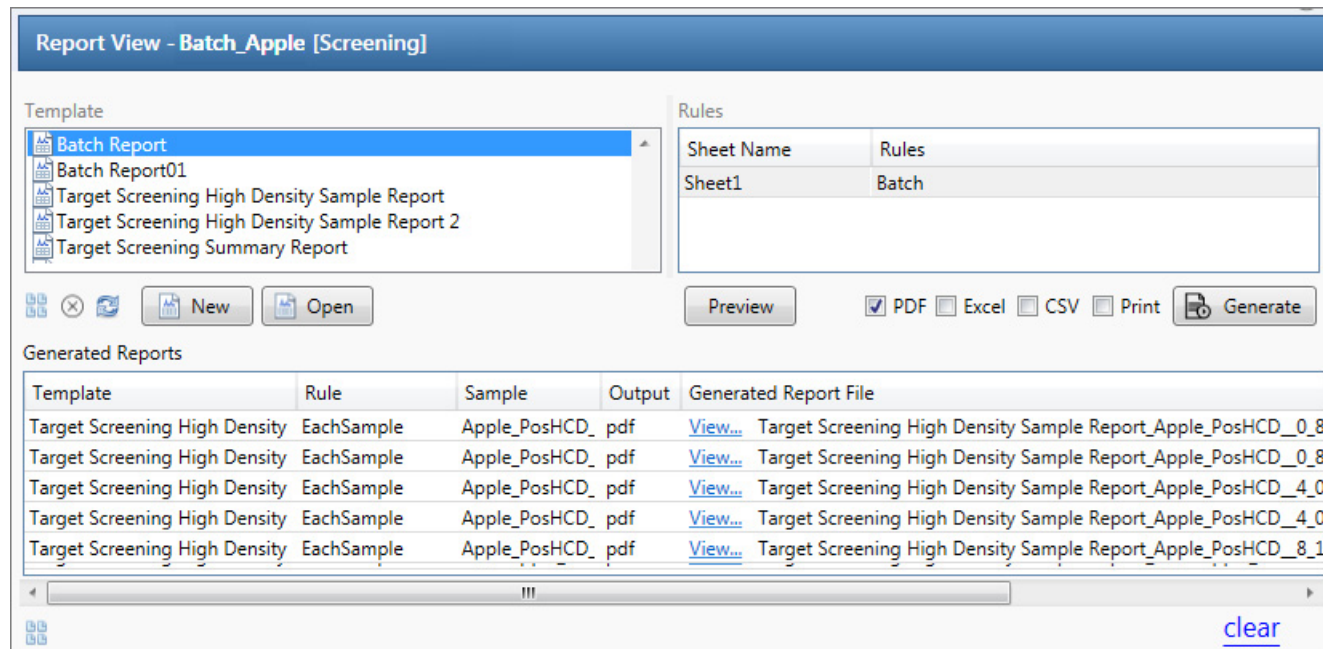
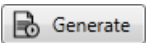



Table 57. Report View parameters (Sheet 1 of 2)

Parameter	Description
Template	
	Displays all report templates.
Rules	
Sheet Name	Specifies each sheet in the report.
Rules	Specifies the type of data used in each sheet in the selected report. <ul style="list-style-type: none">• Batch• EachSample• SampleType: <i>SampleType</i>• CompoundType: <i>CompoundType</i>• SampleCustomFormula:
Buttons	
View Report Templates	Displays the C:\TraceFinderData\4.0\Templates\ReportTemplates folder that contains all report templates.
Open	Opens the selected report template in the Report Designer.
New	Opens a blank report template in the Report Designer.
Preview	Opens the Report Designer showing the report information for the current batch in the selected report template format.

Table 57. Report View parameters (Sheet 2 of 2)

Parameter	Description
PDF	Writes the generated report to a PDF file in the ...\\TraceFinderData\\4.0\\Projects\\ <i>batch</i> \\ReportOutput folder.
Excel	Writes the generated report to a PDF file in the ...\\TraceFinderData\\4.0\\Projects\\ <i>batch</i> \\ReportOutput folder.
CSV	Saves the generated report as a PDF file in the ...\\TraceFinderData\\4.0\\Projects\\ <i>batch</i> \\ReportOutput folder. When the report contains multiple sheets, the application writes each sheet as a separate CSV file.
Print	Prints the generated report to your default printer.
 Generate	Generates the selected type of reports for the current batch using the selected report template.
Generated Reports	
Template	Report template used for the report. See Example of a report template list .
Rule	Type of data used in each sheet of the report. See Rules .
Sample	For sample-level reports, the name of each sample in the report.
Output	Type of output specified for the report: PDF, Excel, CSV, or Print.
Generated Report File	Lists the output file name for each report in the ...\\TraceFinderData\\4.0\\Projects folder.
View	Displays the generated output file.
 View Generated Reports	Displays the C:\\TraceFinderData\\4.0\\Projects folder that contains all report outputs.
Clear	Removes all reports from the Generated Reports display. This does not delete the reports from the C:\\TraceFinderData\\4.0\\Projects folder.

Working in the Local Method View for Target Screening Batches

A local method is a copy of a master method associated with a batch. You can edit only the local copy of the method, or you can edit the master method and overwrite the local copy with the edited master method.

In the Local Method view, you can edit the local method parameters. A local method is a copy of a master method associated with a batch. Local methods are named *Batch_MasterMethod*.

❖ To open the Local Method View

1. Click **Analysis** in the navigation pane.
2. Click **Local Method**.



The [Local Method View](#) for the currently selected batch opens.

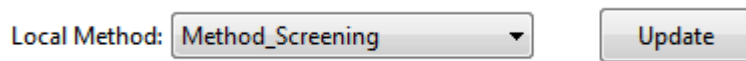
You can edit many of the method parameters in a local method. Editing the local method does not affect parameters in the master method.

For detailed descriptions of method parameters, refer to “Editing a Master Method” in Chapter 5, “Using the Method Development Mode for Target Screening Methods,” in the *TraceFinder Lab Director User Guide*.

3. Enter any local changes to the method.
4. When you have finished editing the local method, choose **File > Save**.
5. To process the batch or create new reports with the edited local method, return to the Batch View and submit the batch.

❖ To overwrite the local method with the master method in the Batch View

In the Batch View, click **Update**.



The application overwrites the local method with the master method of the same name. You can use this feature to overwrite an edited local method with the original master method or to overwrite the local method with an updated master method.

Figure 112. Local Method View

Analysis ▾		Local Method View - Batch_4_10_screening_Method_screening	
▶ Batch View		Master method: Method_screening	
▶ Data Review			
Report View			
▼ Local Method ▶			
Acquisition		Lab Name: <input type="text" value="Default Laboratory"/>	
Target Screening		Assay type: <input type="text" value="Assay name"/>	
Processing		Injection Volume: <input type="text" value="1.00"/> <input type="button" value="▲"/> <input type="button" value="▼"/>	
Peak Detection		Mass Precision: <input type="text" value="2.00"/> <input type="button" value="▲"/> <input type="button" value="▼"/>	
Reports		Instrument Method: <input type="text" value="Instrument1"/> <input type="button" value="Edit"/> <input type="button" value="Update"/>	
		Notes <input type="text"/>	

5 Using the Analysis Mode for Target Screening Batches

Working in the Local Method View for Target Screening Batches

Using the Analysis Mode for Unknown Screening Batches

Use the features of the Analysis mode to do the following:

- Create unknown screening batches.
- Submit unknown screening batches for acquisition, processing, or report generation.
- Review unknown screening batches, batch data, reports, and local methods.

IMPORTANT TraceFinder 4.1 uses the same data as TraceFinder 4.0. By default, the application stores the method, project, template, and compound database data for the 4.1 release in the TraceFinderData\4.0 folder.

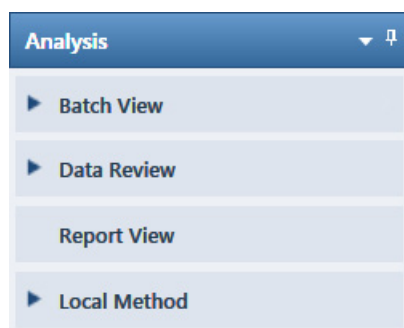
Contents

- [Working in the Batch View for Unknown Screening Batches](#)
- [Working in Data Review for Unknown Screening Batches](#)
- [Working in the Report View for Unknown Screening Batches](#)
- [Working in the Local Method View for Unknown Screening Batches](#)

❖ To access the Analysis mode

Click **Analysis** in the navigation pane.

The Analysis navigation pane opens.



Working in the Batch View for Unknown Screening Batches

In the Batch View, you can manually create and edit a new unknown screening batch or open and edit a previously saved batch. When you submit a batch, you can acquire and process data and optionally create reports for the submitted samples.

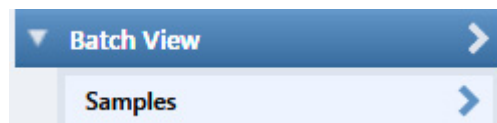
The Analysis mode includes a toolbar:



Use the [Toolbar](#) or the equivalent commands in the [Batch View Shortcut Menu](#) to create the sample list and submit samples for acquisition.

❖ To open the Samples page in the Batch View

1. Click **Analysis** in the navigation pane of the current mode.
2. Click the **Batch View** down arrow and then click **Samples**.



The Samples page opens.

Related Topics

- [Samples Page Features](#)
- [Creating a New Batch](#)
- [Editing a Batch](#)
- [Submitting a Batch](#)

Samples Page Features

Use the Samples page to create a batch.

Batch View - Batch_A														
Local Method:		Method_Unknown	Update		Instrument:		Thermo Scientific Instrument		User:					
	Status	Filename	Sample type	Groups	Level	Sample ID	Sample name	Vial position	Injection volume	Conversion Factor	Channel	Barcode Expected	Barcode Actual	Sample Volume
▶ 1	●	Unknown1	Specimen						10.0	1.000	Auto			1
2	●	Unknown2	Specimen						10.0	1.000	Auto			1
3	●	Unknown3	Specimen						10.0	1.000	Auto			1

Samples Page

The Samples page includes the following features:

- [Column Display](#)
- [Status Indicators](#)
- [Groups](#)
- [Sample Weight Calculation](#)
- [Instrument Methods](#)
- [Toolbar](#)
- [Batch View Sample List](#)
- [Batch View Shortcut Menu](#)

Column Display

The sample list contains many columns of information. You can scroll to see all the columns, and you can customize which ones to display and their display order.

❖ To scroll the sample list

Use the horizontal scroll bar at the bottom of the sample list to view all the information.

When you use the scroll bar at the bottom of the sample list, the following columns remain fixed: Status, Filename, Sample Type, Groups, Level, Sample ID, and Sample Name, while the other columns scroll right and left.

❖ To customize the column display

1. Right-click the sample list and choose **Modify Columns**.

The [Modify Columns dialog box](#) opens.

2. Use the arrow buttons to move all the columns that you want displayed to the Displayed Columns pane.

These columns appear after the Status, Filename, Sample Type, Groups, Level, Sample ID, and Sample Name columns.

3. To arrange the order of the columns, do the following:

- a. In the Displayed Columns pane, select a column name.
- b. Use **Up** or **Down** to move the selected column up or down in the list.

The first column in the list represents the leftmost column in the Batch View sample list, and the last column in the list represents the rightmost column in the Batch View sample list.

Note The following columns are fixed: Status, Filename, Sample Type, Groups, Level, Sample ID, and Sample Name.

6 Using the Analysis Mode for Unknown Screening Batches

Working in the Batch View for Unknown Screening Batches

4. To change the width of a column, do the following:
 - a. In the Displayed Columns pane, select the column width.

5	Sample ID	100
▶ 6	Sample name	100
7	Vial position	100

- b. Type a new value for the width.
5. Repeat [step 4](#) for all columns whose widths you want to change, and click **OK**.

The columns in the sample list immediately reflect your changes. The application uses these settings for all sample lists in the Batch View.

Figure 113. Modify Columns dialog box

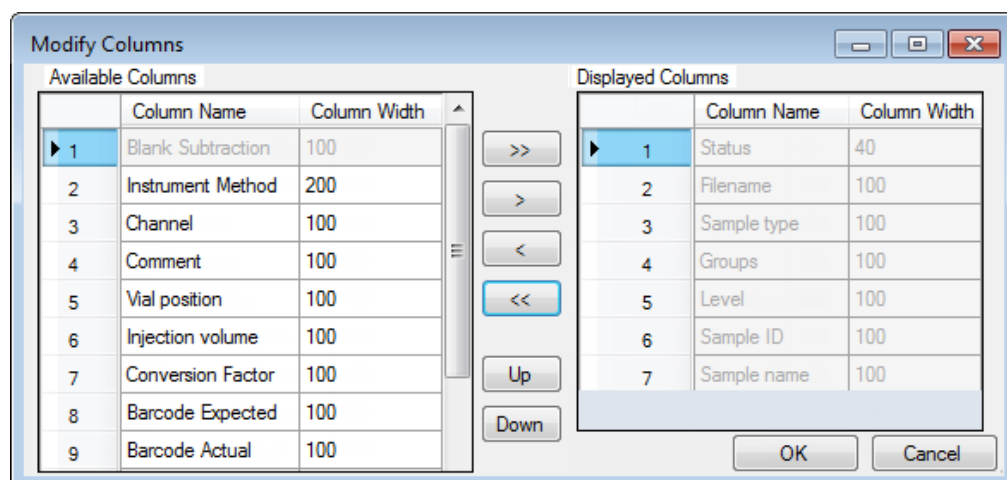






Table 58. Button descriptions for the Modify Columns dialog box




Button	Description
	Moves all columns to the Displayed Columns pane.
	Moves the selected column to the Displayed Columns pane.
The following buttons apply to all columns, except for those that are fixed: Status, Filename, Sample Type, Groups, Level, Sample ID, and Sample Name.	
	Moves the selected column to the Available Columns pane.
	Moves all columns except fixed columns.
	Moves the selected column name in the Displayed Columns pane one row up in the column order.
	Moves the selected column name in the Displayed Columns pane one row down in the column order.

Status Indicators

Status indicators show the current status of each sample during the acquisition and processing.

-  Sample is not acquired.
-  Sample is acquired but not processed.
-  Sample is acquired and processed.
-  Sample is currently acquiring.

Local Method: Method_Alprazolam Update

	Status	Filename	Sample type	Groups	Sample ID	Sample name	Vp
1		Solvent_A	Solvent				1
2		Cal_5	Calibrator				2
3		Unknown1	Specimen				3

Status indicators




Groups

Use the Groups feature to assign samples to a group.

❖ To create a group

- For each sample, type the name of a group in the Groups column.
Repeat this for each sample that you want to include in a group.
- Create as many groups as you want.

Note To assign a sample to multiple groups, separate the groups with a comma.

Status	Filename	Groups
	Benzo26473	groupB, groupA
	Benzo25557	groupB
	Benzo26154	groupB, groupA

The application uses the first group listed for the Maximum Fold analysis and in the Group Averages pane.

You can create a control group (specify the group name as Control) to use for group averages in Data Review.

For information about group averages, see [Group Averages Pane](#). For information about maximum fold analysis, see [Cross Sample Peak List Pane](#).

Sample Weight Calculation

Use the sample weight features to calculate the conversion factor for a sample. The application uses different methods to calculate the conversion factor for liquid or solid calculation types.

Liquid: $SampleVolume \div DilutionFactor$

Solid: $(SampleVolume \times DilutionFactor) \div SampleWeight$

Manual: The application does not calculate the Conversion Factor. Instead, you can enter the Conversion Factor value.

Follow these procedures:

- To display the features for calculating sample weight
- To calculate the conversion factor for a liquid sample
- To calculate the conversion factor for a solid sample
- To manually specify the conversion factor for a sample

❖ To display the features for calculating sample weight

If the Conversion Factor, Sample Volume, Dilution Factor, Sample Weight, Calculation Type, and Final Units columns are not visible, right-click and choose **Enable Sample Weight Calculation**.

Conversion Factor	Sample Volume	Dilution Factor	Sample Weight	Calculation Type	Final Units
1.000	1	1	1	Liquid ▼	
1.000	1	1	1	Solid ▼	
1.000	1	1	1	Manual ▼	

❖ To calculate the conversion factor for a liquid sample

Note The application uses the following formula to calculate the Conversion Factor:
 $SampleVolume \div DilutionFactor$

1. From the **Calculation Type** list, select **Liquid**.
For a liquid sample, the Sample Weight value is not editable.
2. In the Sample Volume column, type the volume in ng/mL for your sample.
3. In the Dilution Factor column, type the value for the dilution.

For example, if you have 1000 ng/mL of a substance that is too concentrated for the mass spectrometer, you can dilute it by 1000. Then your injection volume is 1, your conversion factor is 1000, and your sample amount is 1000.

4. In the Final Units column, type the units that you want to use for the calculated amount in the Data Review view or in reports.

❖ **To calculate the conversion factor for a solid sample**

Note The application uses the following formula to calculate the Conversion Factor:
 $(SampleVolume \times DilutionFactor) \div SampleWeight$

1. From the **Calculation Type** list, select **Solid**.
2. In the Sample Weight column, type the weight in ng for your sample.
3. In the Sample Volume column, type the volume in ng/ml for your sample.
4. In the Dilution Factor column, type the value for the dilution.

For example, if you have 1000 ng/ml of a substance that is too concentrated for the mass spectrometer, you can dilute it by 1000. Then your injection volume is 1, your conversion factor is 1000, and your sample amount is 1000.

5. In the Final Units column, type the units that you want to use for the calculated amount in the Data Review view or in reports.

❖ **To manually specify the conversion factor for a sample**

Note The application uses the specified conversion factor when it calculates the amount for the sample.


1. From the **Calculation Type** list, select **Manual**.
 For a manually calculated sample, the only available columns are the Conversion Factor and the Final Units.
2. In the Conversion Factor column, type the conversion factor to use for your sample.
3. In the Final Units column, type the units that you want to use for the calculated amount in the Data Review view or in reports.

Instrument Methods

Use the Instrument Methods column to specify instrument methods for the samples.

Note By default, the Instrument Method column is not displayed in the Batch View sample list.

❖ **To specify instrument methods for samples**

1. Display the Instrument Method column in the sample list:
 - a. Right-click the sample list and choose **Modify Columns**.
 The Modify Columns dialog box opens.
 - b. In the Available Columns pane, select **Instrument Method**.
 - c. Click  to move the Instrument Method column to the Displayed Columns pane.

6 Using the Analysis Mode for Unknown Screening Batches

Working in the Batch View for Unknown Screening Batches

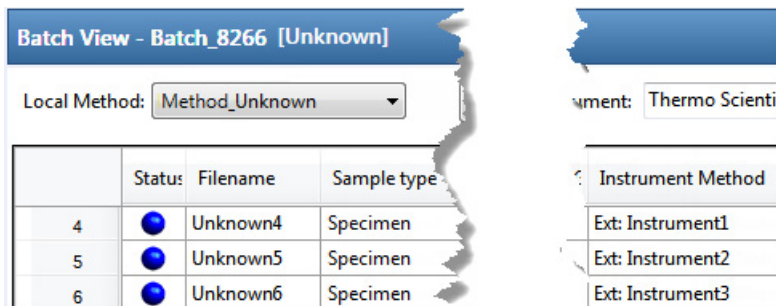
d. Click **OK**.

The application displays the Instrument Method column, defaulting to the instrument method specified in the master method.

2. Click the Instrument Method column and select an instrument method from the list.

This list contains all the available instrument methods. Instrument methods from external sources are prefixed with “Ext:”.

You can specify a different instrument method for each sample.



When you submit the batch for acquisition, the application saves a copy of the selected instrument methods to the following folders:

External instrument methods:

...\\TraceFinderData\\4.0\\Projects\\...\\batch\\Methods\\method\\ExternalMethods

Local instrument methods:

...\\TraceFinderData\\4.0\\Projects\\...\\batch\\Methods\\method

Toolbar








The Analysis mode includes this toolbar for creating and submitting a batch.



Table 59. Toolbar icons (Sheet 1 of 2)

Icon	Description
	Adds the specified number of new, empty samples to the end of the sample list. See the instructions To add samples to the list .
	Inserts a new, empty sample or samples above the selected sample. See the instructions To insert samples into the list .
	Removes the selected samples from the sample list. See the instructions To remove samples from the list .
	Adds imported samples from a CSV, an XML, or an SLD file to the sample list. See the instructions To import samples into the list .

Table 59. Toolbar icons (Sheet 2 of 2)

Icon	Description
	Submits only the selected samples for acquisition, processing, or report generation. See the instructions To submit samples in the batch.
	Submits the batch for acquisition, processing, or report generation. See the instructions To submit samples in the batch.
	Submits only the selected samples for processing. See the instructions To submit selected peaks for processing.
	Opens the Acquisition mode where you can use a batch template to define a standard sequence composed of various sample types to be assembled into a batch of samples. See Working in Data Review for Unknown Screening Batches.
	Opens the Acquisition mode where you can create a batch template that contains the basic settings and sample types for your batches. See Using the Acquisition Mode.
	Opens the Quick Acquisition window where you can quickly submit a single sample. See Appendix A, "Using Quick Acquisition."
	Opens the Audit Viewer where you can view audit logs. See Chapter 8, "Using the Audit Viewer." Available only when you enable Auditing in the Administrator Console. Refer to the instructions in the <i>TraceFinder Administrator Console User Guide</i> .

Batch View Sample List

The sample list displays all the quantitative data for the samples of a batch. Status indicators for each sample indicate if the sample is currently acquiring, not acquired, acquired, or processed.

The sample list includes the following columns of information:

Figure 114. Batch View sample list




Local Method:

Method_Unknown

Update

Instrument:

Thermo Scientific Instrument

	Status	Filename	Sample type	Groups	Level	Sample ID	Sample name	Vial position	Injection volume
1		UnknownA1	Specimen			1		CStk1-01:7	10.0
2		UnknownA2	Specimen			1		CStk1-01:7	10.0
3		UnknownA3	Specimen			1		CStk1-01:7	10.0

Calculation Type	Conversion Factor	Dilution Factor	Sample Weight	Sample Volume	Final Units
Liquid <div></div>	1.000	1	1	1	
Liquid <div></div>	1.000	1	1	1	
Liquid <div></div>	1.000	1	1	1	

Instrument Method	Channel	Barcode Expected	Barcode Actual	Comment
Instrument1 <div></div>	Auto			
Instrument1 <div></div>	Auto			
Instrument1 <div></div>	Auto			

Table 60. Batch View sample list columns (Sheet 1 of 3)







Column	Description
Status	 Sample is not acquired.  Sample is acquired but not processed.  Sample is acquired and processed.  Sample is currently acquiring.
Note When you include unknown screening features in a quantitation method and you choose to process with only the unknown screening criteria, both the Sample View and the Unknown Screening View show the Status for the samples as acquired and processed ().	
Note When you include unknown screening features in a target screening method and you choose to process with only the unknown screening criteria, both the Samples view and the Unknown Screening View show the Status for the samples as acquired and processed ().	

Table 60. Batch View sample list columns (Sheet 2 of 3)

Column	Description
Filename	Name of the raw data file that contains the sample data.
Sample Type	Defines how the application processes the sample data. Each sample is classified as one of the following sample types: Specimen, QC, Solvent, Hydrolysis, Unextracted, or Negative. Default: Specimen
	Note Unknown Screening does not use calibration sample types (Calibrator).
Groups	Threshold group to which a sample belongs.
Level	Not used in Unknown Screening processing.
Sample ID	A user-defined, alphanumeric string that identifies a sample.
Sample Name	A user-defined name that identifies a sample.
Vial Position	The tray vial number used for an autosampler acquisition.
Injection Volume	The injection volume (in microliters) of the injected sample. When you are using an autosampler, you can set the default injection volume in the Autosampler dialog box in the Instrument View. The minimum and maximum injection volumes that you can use depend on the Autosampler you configure. The usable range depends on the injection mode and might be smaller than the displayed range. The Injection Volume value set in the master method overwrites the value in the instrument method. Valid range: 0.1 through 5000 µL
Calculation Type	Liquid: The application calculates the Conversion Factor as $\text{SampleVolume} \div \text{DilutionFactor}$ Solid: The application calculates the Conversion Factor as $(\text{SampleVolume} \times \text{DilutionFactor}) \div \text{SampleWeight}$ Manual: Sample Volume, Dilution Factor, Sample Weight, and Final Units columns are not available, and the Conversion Factor value is editable.
Conversion Factor	Editable only when Calculation Type is Manual. Default: 1
Sample Volume	Default: 1
Dilution Factor	Default: 1
Sample Weight	Available only when Calculation Type is Solid. Default: 1

Table 60. Batch View sample list columns (Sheet 3 of 3)

Column	Description
Final Units	Specifies the calculated amount in the Data Review view or in reports. Default: 1
Instrument Method	Specifies the instrument to use for the acquisition. This column is hidden by default. To display this column, see To customize the column display .
Channel	Specifies the channel on which the sample was run. If the sample is not acquired, the value is Pending. The Channel column is available only when you have activated multiplexing in the Configuration console. Refer to “Multiplexing” in Chapter 1, “Using the Configuration Console,” in the <i>TraceFinder Lab Director User Guide</i> .
Barcode Expected	A user-entered barcode for the vial.
Barcode Actual	An actual barcode for the vial. This value is not editable.
Comment	A user-defined comment for the sample.

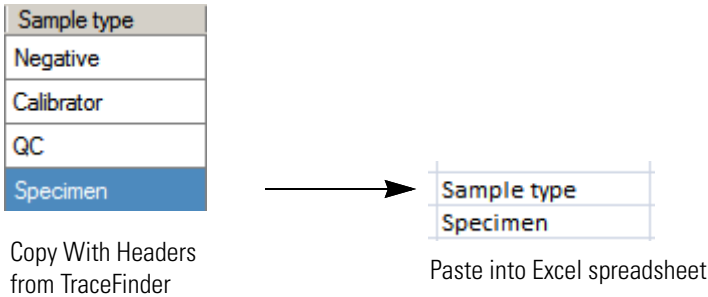
Batch View Shortcut Menu

The Batch View includes a shortcut menu for creating a batch.

Table 61. Batch View shortcut menu commands (Sheet 1 of 2)

Command	Description
Add Sample	Adds a single empty row to the sample grid.
Insert Sample	Inserts a single empty row to the sample grid above the selected row.
Insert Copy Sample	Copies the currently selected row and inserts a copy above the row.
Reinject Selected Samples	Creates a copy of the selected sample and appends INJ001 to the file name. Additional reinjections of the same sample are numbered INJ002, INJ003, and so forth.
Remove Selected Samples	Removes selected samples from the sample grid.
Import Samples	Opens the Sample Import Tool. See To import samples into the list .
Browse in Raw File (Move)	Opens a dialog box where you can select a raw data file to use for the selected sample row. The application removes the raw data file from the source location.
Browse in Raw File (Copy)	Opens a dialog box where you can select a raw data file to use for the selected sample row. The application copies the raw data file from the source location.
Map Raw Files to Samples	Opens a dialog box where you can select multiple raw data files to use for the selected sample rows.
Copy Down	Copies the value in the selected row to all rows below it. This command is available only when you have selected a value that can be copied down.

Table 61. Batch View shortcut menu commands (Sheet 2 of 2)

Command	Description
Fill Down	Enters sequential values in the column starting with the value in the selected row and ending with the last row in the column. This command is available only when you have selected a value that can be filled down.
Modify Columns	Opens the Modify Columns dialog box .
Enable/Disable Sample Weight Calculation	Displays or hides the Sample Volume, Dilution Factor, Sample Weight, Calculation Type, and Final Units columns.
Copy	Copies the data in the selected rows or columns to the Clipboard. Use this command to copy sample information into a text editor or spreadsheet application. You cannot paste this data back into the Batch View sample list.
Copy with Headers	Copies the data in the selected rows or columns and the associated column headers to the Clipboard. Use this command to copy sample information into another text editor or spreadsheet application. You cannot paste this data back into the sample list.
<p>For example</p>  <p>Copy With Headers from TraceFinder</p> <p>Paste into Excel spreadsheet</p>	
Paste	Pastes a single column of copied data from another text editor or spreadsheet application into the selected column.
Undo Last Paste	Removes the last pasted item in the Batch View.
Export to CSV File	Opens the Save As dialog box where you can save the current sample list to a CSV file.
Edit Instrument Method	<p>Opens the Instrument Setup window where you can edit the parameters of the instrument method.</p> <ul style="list-style-type: none"> When you edit an external method, the application updates the method in the ...\\Xcalibur\\methods folder. When you edit an internal method, the application updates the method in the ...\\TraceFinderData\\4.0\\Projects\\...\\batch\\Methods\\method folder. <p>For detailed information about editing instrument methods, refer to Chapter 3, “Using Instrument Methods in the Method Development Mode,” in the <i>TraceFinder Lab Director User Guide</i>.</p>

Creating a New Batch

In the Batch View, you can create a new batch.

Follow these procedures:

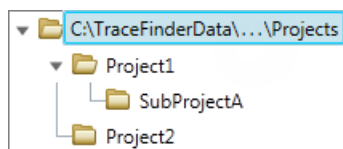
- [To create a new batch](#)
- [To add samples to the list](#)
- [To insert samples into the list](#)
- [To import samples into the list](#)
- [To remove samples from the list](#)
- [To copy a sample](#)
- [To reinject a sample](#)
- [To edit sample values](#)
- [To browse in raw data files](#)
- [To customize the column display](#)

❖ To create a new batch

1. Choose **File > New > Batch** from the main menu.

The [Create New Batch Dialog Box](#) opens, displaying all drives that contain projects.

2. Select a drive from the list.



Tip The application displays all configured and enabled repositories.

3. Select the folder where you want to store your batch.

Tip To activate the Create button, you must enter a unique batch name. If the Create button is not activated, you have entered a batch name that is already used.

To create a new folder for the storage location, see [Editing Folders for Batches](#).

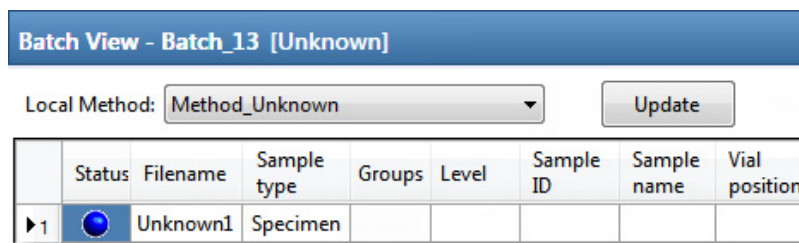
4. Select **Unknown Only** from the Type list.

The batch list displays all batches in the selected folder. The Method list displays all methods for the unknown screening type.

5. Select a master method from the Master Method list.

6. Click **Create**.

A new batch opens with one Specimen sample.



The batch name in the title bar indicates that you are creating an unknown screening batch.

❖ **To add samples to the list**

1. To add a single sample row, right-click the sample list and choose **Add Sample**.
2. To add multiple sample rows, select the number of rows and then click the **Add Sample** icon, .

The application adds the specified number of new, empty samples to the end of the sample list.

❖ **To insert samples into the list**

Select the sample above which you will insert new, Specimen samples, and then do one of the following:


- To insert a single sample row, right-click and choose **Insert Sample**.
- To insert multiple sample rows, select the number of rows and then click the **Insert Sample** icon .

The application inserts the Specimen samples above the selected sample.

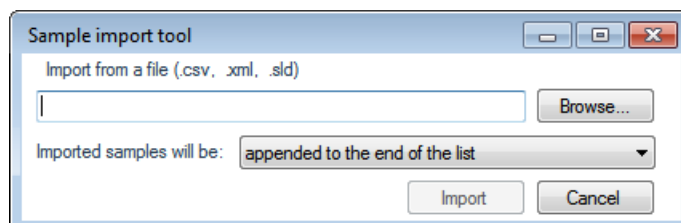
Inserted samples

	Status	Filename	Sample type	Groups	Level
1		cal_std_5	Calibrator		
2		Unknown2	Specimen		
3		Unknown1	Specimen		
4		cal_std_10	Calibrator		

❖ To import samples into the list

1. Choose **Batch > Import Samples** from the main menu, or click the **Import Samples** icon, .

The Sample Import Tool dialog box opens.



From this dialog box, you can import samples from a CSV, an XML, or an SLD file.

2. Click **Browse** and select a CSV, an XML, or an SLD file that contains the samples to import.
3. From the Imported Samples Will Be list, select **Appended to the End of the List** or **Inserted at the Selected Row**.
4. Click **Import**.

The Sample Import Tool dialog box closes, and the application adds the specified samples to the sample list.

When you import samples from an Xcalibur sequence file (.sld), the application makes the following column name substitutions:

Xcalibur column	TraceFinder column
Position	Vial Position
Inj Vol	Injection Volume
Dil Factor	Conversion Factor

When you import samples from an Xcalibur sequence file (.sld), the application makes the following sample type substitutions:

Xcalibur sample type	TraceFinder sample type
Blank	Negative
Std Bracket	Calibrator

5. (Optional) When using multiplexing, select a channel for each imported sample. Imported samples default to Auto.

Note The Channel column is available only when you have activated multiplexing in the Configuration console. Refer to “Multiplexing” in Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.

❖ **To remove samples from the list**

1. Select the samples that you want to remove.

Tip Use the CTRL or SHIFT keys to select multiple samples.

2. Right-click and choose **Remove Selected Samples**.

❖ **To copy a sample**

1. Select the sample that you want to copy.
2. Right-click and choose **Insert Copy Sample**.

The application inserts the copy above the selected sample.

❖ **To reinject a sample**

1. In the sample list, select the sample that you want to reinject.
2. Right-click and choose **Reinject This Sample**.

The application creates a copy of the selected sample and appends INJ001 to the file name. Additional reinjections of the sample are numbered INJ002, INJ003, and so forth. The application copies all parameter values from the original sample.

❖ **To edit sample values**

1. For each sample, do one of the following:

Type a new file name over the current filename.

—or—

Double-click the Filename column and locate a raw data file to use for the sample.

—or—

Right-click and choose **Browse in Raw File**, and then locate a raw data file to use for the sample.

By default, the application sets the Sample Type to Unknown.

2. For each sample, click the Sample Type column and select a sample type from the list.

Available sample types			
Specimen	Hydrolysis	Solvent	QC
Unextracted	Negative		

3. Type a vial position in the Vial Position column for each sample.
4. Type a volume in the Injection Volume column for each sample.

The minimum injection volume value allowed is 0.1 µL; the maximum injection volume value allowed is 5000 µL.

5. (Optional) Type or edit the values for the remaining columns.

Note When you use the scroll bar at the bottom of the sample list, the following columns remain fixed: Status, Filename, Sample Type, Groups, Level, Sample ID, and Sample Name, while the other columns scroll right and left.

To automatically copy or fill values in these columns, see [Appendix B, “Using Copy Down and Fill Down.”](#)

❖ To browse in raw data files

1. Do one of the following:

Double-click the Filename column.


–or–

Right-click and choose **Browse in Raw File**.

The What Raw File Would You Like to Use dialog box opens.

2. Select a raw data file to use for the sample or use the CTRL key to select multiple files, and then click **Open**.

The application overwrites the selected, unacquired sample in the batch with the first “browsed in” file and adds any additional browsed in files below the selected sample.

For all browsed-in raw data files, the application sets the Status to Acquired, , and sets the Sample Type to Specimen.

Note You cannot overwrite an acquired sample. When you select a sample that is acquired, the application adds all browsed in files below the selected sample.

❖ To customize the column display

1. Right-click the Batch View sample list and choose **Modify Columns**.

The [Modify Columns dialog box](#) opens.

2. Use the arrow buttons to move all the columns that you want displayed to the Displayed Columns pane.

All the columns you select appear after the Status, Filename, Sample Type, Groups, Level, Sample ID, and Sample Name columns.

3. To arrange the order of the columns, do the following:

- a. In the Displayed Columns pane, select a column name.

- b. Use **Up** or **Down** to move the selected column up or down in the list.

The first column in the list represents the leftmost column in the Batch View sample list, and the last column in the list represents the rightmost column in the Batch View sample list.

Note The following columns are fixed: Status, Filename, Sample Type, Groups, Level, Sample ID, and Sample Name.

4. To change the width of a column, do the following:
 - a. In the Displayed Columns pane, select the column width.

7	Vial position	69
8	Injection volume	100
9	Calculation Type	100

- b. Type a new value for the width.
5. When you have completed your changes, click **OK**.

The columns in the sample list immediately reflect your changes. The application uses these settings for all sample lists in the Batch View.

Figure 115. Modify Columns dialog box

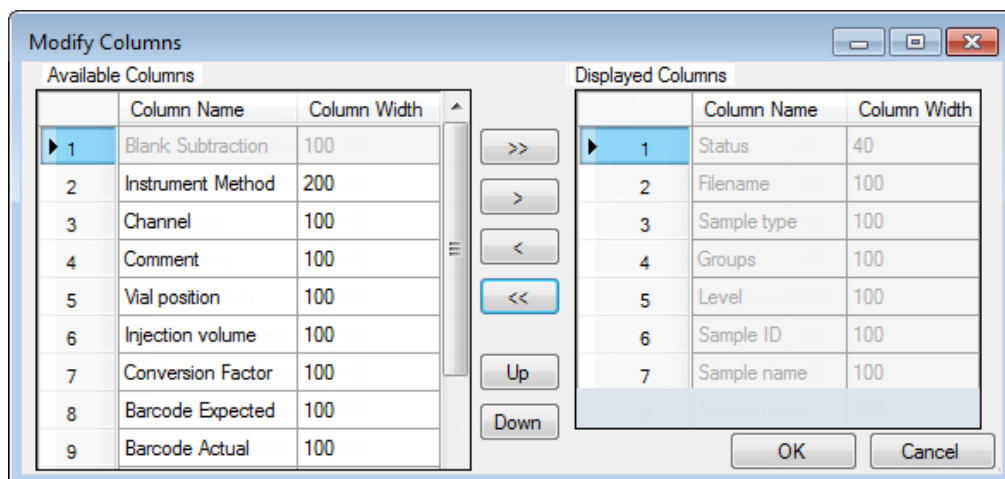


Table 62. Button descriptions for the Modify Columns dialog box

Button	Description
	Moves all columns to the Displayed Columns pane.
	Moves the selected column to the Displayed Columns pane.
The following buttons apply to all columns, except for those that are fixed: Status, Filename, Sample Type, Groups, Level, Sample ID, and Sample Name.	
	Moves the selected column to the Available Columns pane.
	Moves all columns except those that are fixed.
	Moves the selected column name in the Displayed Columns pane one row up in the column order.
	Moves the selected column name in the Displayed Columns pane one row down in the column order.

Create New Batch Dialog Box

Use the Create New Batch dialog box to select a folder and method for your batch and to name the new batch.

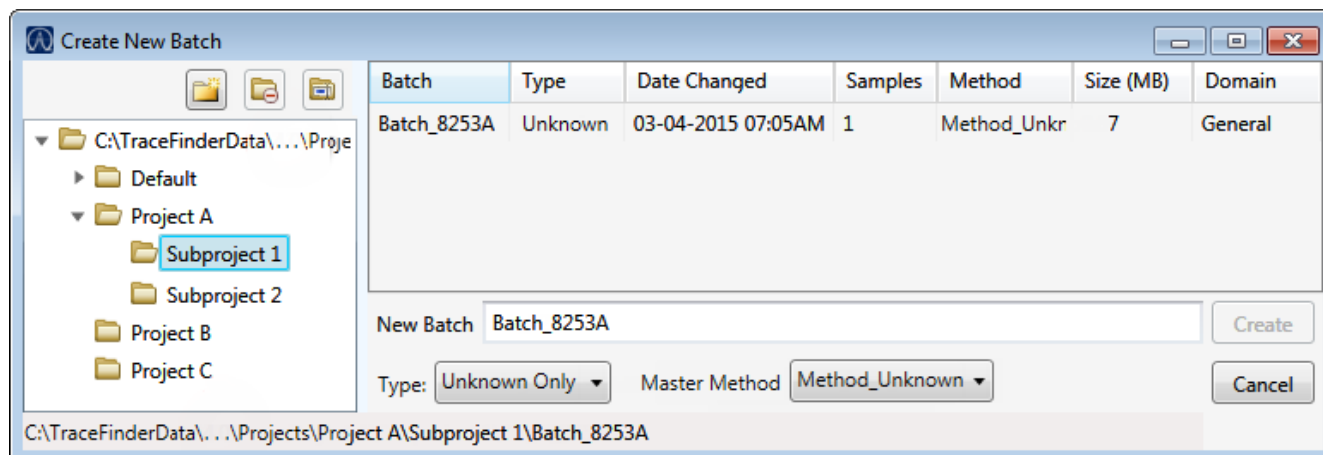


Table 63. Create New Batch dialog box parameters (Sheet 1 of 2)




Parameter	Description
 Create New Folder	Adds one of the following: <ul style="list-style-type: none"> When a drive is selected, adds a new project-level folder to the drive. When a project folder is selected, adds a subproject-level folder to the selected project. When a subproject folder is selected, adds a lower-level folder to the subproject. Or, you can right-click and choose Create Folder.
 Delete Folder	With no confirmation prompt, immediately removes the selected folder. You cannot delete a folder that contains lower-level folders; you must delete the lower-level folders first. Or, you can right-click and choose Delete.
 Rename Folder	Renames the selected folder. Or, you can right-click and choose Rename.
Batch table	
Batch	Name of batches in the selected project.
Type	Type of batch: Quan, Screening, or Unknown Only.
Date Changed	Date that the batch was last updated.
Samples	Number of samples in the batch.
Method	Name of the method used to create the batch.
Size	Size of the batch in megabytes.
Domain	TraceFinder domain in which the batch was created.

Table 63. Create New Batch dialog box parameters (Sheet 2 of 2)

Parameter	Description
New batch parameters	
New Batch	Name of the new batch to create. Note If the Create button is not activated, you have entered a name that is already used or you have not selected a method.
Type	Type of batch to create: Quan, Screening, or Unknown Only.
Method	Method used to create the new batch.
Path	Path to the project in the TraceFinderData\4.0\Projects folder where the batch is created.
Buttons	
Create	Creates the specified batch and opens the Batch View for the new batch.
Cancel	Closes the Create New Batch dialog box without creating a batch.


Editing Folders for Batches

From the Create New Batch dialog box, you can create new folders for your batches. You can also delete or rename folders.

Use these procedures:

- [To create new project folders](#)
- [To delete project folders](#)
- [To rename project folders](#)

❖ To create new project folders

1. In the Create New Batch dialog box, select the folder for which you will create a new lower-level folder.
 - You can select the main TraceFinderData\4.0\Projects folder and create a new folder under it.
 - You can select one of the existing folders and create a lower-level folder under it.
2. Click the **Create Folder** icon, .


The application adds a new lower-level folder to the selected folder.

3. Select the new folder name and type a name for the folder.

Folder names are limited to 30 characters and can contain spaces and special characters, except for the following special characters: \ / : + ? " < >

Note After you add a lower-level folder, you cannot rename the parent folder.


❖ To delete project folders

1. In the Create New Batch dialog box, select the folder to delete.
2. Click the **Delete Folder** icon, .

With no confirmation prompt, the application immediately removes the selected folder.

Note You cannot delete folders that contains lower-level folders; you must delete the lower-level folders first.

❖ To rename project folders

1. In the Create New Batch dialog box, select the folder to rename.
2. Click the **Rename Folder** icon, .

Note You cannot rename folders that contain lower-level folders.

3. Type a new name for the folder and press ENTER.

The application saves the new folder name.

Editing a Batch

In the Batch View, you can open a saved batch and edit the sample list. You can add samples, edit samples, or remove samples. If the batch has already been acquired, you can select specific samples for reinjection. If the batch has unacquired samples when you complete your edits, you can save it as a “ready to acquire” batch.

Follow these procedures:

- [To open a saved batch](#)
- [To open a recent batch](#)
- [To edit samples in a batch](#)
- [To reinject a sample from a previously acquired batch](#)

❖ To open a saved batch

1. Choose **File > Open > Batch** from the main menu.

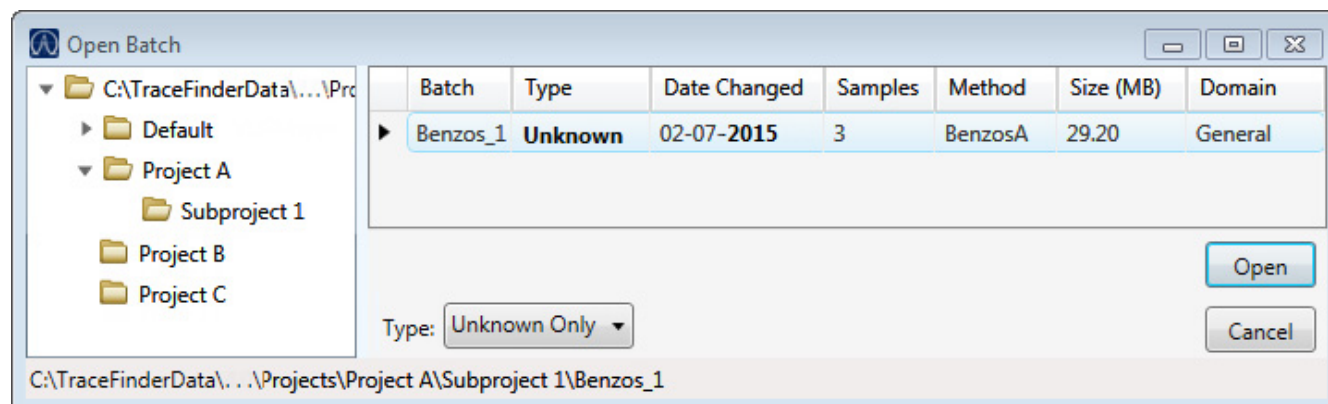
The [Open Batch dialog box](#) opens.

2. Select a project and a subproject.
3. Select **Unknown Only** or **Any** from the Type list.

The batch list displays all batches created with Unknown methods (or all methods of all types when you select Any).

4. Select a batch from the list.
5. Click **Open**.

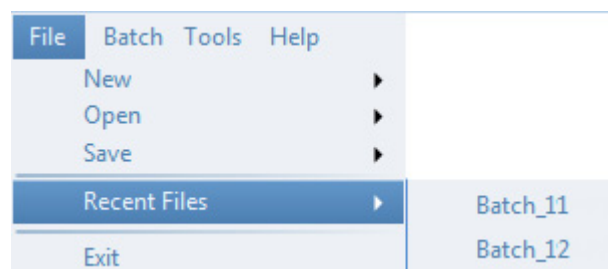
The selected batch opens in the Batch View.

Figure 116. Open Batch dialog box**Table 64.** Open Batch dialog box parameters

Parameter	Description
Batch	Name of batches in the selected project.
Type	Type of batch: Quan, Screening, or Unknown Only.
Date Changed	Date the batch was last updated.
Samples	Number of samples in the batch.
Method	Name of the method used to create the batch.
Size	Size of the batch in megabytes.
Domain	TraceFinder domain in which the batch was created.
Path	Path to the project in the TraceFinderData\4.0\Projects folder where the batch is stored.
Buttons	
Type	Type of batch to display in the Batch list: Quan, Screening, Unknown Only, or Any.
Open	Opens the Batch View for the selected batch.
Cancel	Closes the Open Batch dialog box without opening a batch.

❖ To open a recent batch

Choose **File > Recent Files > *batch*** from the main menu.



The selected batch opens in the Batch View.

❖ To edit samples in a batch

Use the commands described in [Working in the Batch View for Unknown Screening Batches](#).





You can add new samples, edit samples, or delete samples.

❖ To reinject a sample from a previously acquired batch

1. In the sample list, select the sample that you want to reinject.
2. Right-click and choose **Reinject This Sample**.

The application creates a copy of the selected sample and appends INJ001 to the file name. Additional reinjections of the sample are numbered INJ002, INJ003, and so forth. The application copies all parameter values from the original sample.

A green status icon indicates previously acquired samples (acquired and processed), and the sample name is grayed out. A blue status icon indicates samples created for reinjection (not acquired).

	cal_50_INJ001	Calibrator	10
	cal_50	Calibrator	10
	cal_10_INJ001	Calibrator	10
	cal_10	Calibrator	10

When you submit all samples in this batch, the application acquires all samples (including previously acquired samples).

3. To save this batch with the new samples for reinjection, choose **File > Save > Batch** from the main menu.

The batch is saved as a prepared batch that is ready to submit. You can open this batch from the Reinject Samples page in the Acquisition mode and submit the batch. The application acquires only the samples that have not been previously acquired.

Submitting a Batch


In the Batch View, you can submit an entire batch or only selected samples in the batch. When you submit a batch for acquisition and processing, you can choose to create reports for the submitted samples. See [Submit Options dialog box](#).

For a description of commands in the shortcut menu, see [Batch View shortcut menu commands](#).

Follow these procedures:

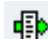

- [To submit selected peaks for processing](#)
- [To submit samples in the batch](#)
- [To view the output files](#)

❖ **To submit selected peaks for processing**

1. In the Peak List, select the **Selected** check box for the peaks that you want to process.
2. Click the **Submit ... for Processing** icon, .

The application processes the selected peaks and updates the data in the Peak Identifications pane.

❖ **To submit samples in the batch**

1. Do one of the following:
 - To submit all samples in the batch, click the **Submit Batch** icon, .
 - To submit specific samples, select the samples and click the **Submit Selected Samples** icon, .

The [Submit Options dialog box](#) opens.

Note You can also submit only selected peaks for processing. See [To submit selected peaks for processing](#).

2. To acquire (or reacquire) the submitted samples, select the **Acquire Data** check box.
 - When all submitted samples have been previously acquired, this option is (by default) not selected.
 - When one or more samples in the batch have not been acquired, this option is (by default) selected.

Tip You can process the data with or without performing peak detection. For example, you might want to turn off peak detection when reprocessing samples.

3. To process the submitted samples, select the **Process Data** check box.
The application displays processing options for the Unknown Screening method.

4. Select the check box for the unknown screening options that you want to use.

Peak Detect: Performs peak detection. You can process the data with or without performing peak detection. For example, you might want to turn off peak detection when reprocessing samples.

Identify: Performs identification. This produces the peak identification data in the Unknown Screening View.

With RT Alignment: Performs retention time alignment. This produces the heat map and group averages data in the Unknown Screening View. When you select this option, the application automatically selects the Peak Detect option.

5. (Optional) Select the **Create Reports** check box.

6. (Optional with multiplexing activated) Select the **Priority Sequence** check box.

The application acquires the priority batch on the next available channel or the assigned channel.

7. (Optional without multiplexing activated) Select the **Priority Sequence** check box and then select one of the following priority options to place the batch in the queue:

- **Next Available Batch** places the batch immediately after the currently acquiring batch.
- **Next Available Sample** places the batch immediately after the currently acquiring sample.

Note When you select Full Sequence Submission in the Configuration console, these options are unavailable because the current batch and the current sample are, in effect, the same thing.

8. To specify the following optional parameters, click **Show Details**.

- a. Select the **Use** check box for the device that you want to use for this acquisition.
- b. Select the **Start Device** check box to indicate the device that will initiate communication with the other instruments.

This is usually the autosampler.

- c. Select the **Start When Ready** check box, which starts all instruments together when they are all ready.

When this is cleared, individual instruments can start at different times and then have to wait for the last instrument to be ready.

- d. Select the system state after it acquires the last batch: **On**, **Standby**, or **Off**.

9. To start the selected processes, click **OK**.

The selected processes begin, and the application shows the real-time display at the bottom of the current window. You can begin another batch in the Analysis mode while you watch the real-time display of the currently acquiring batch.

IMPORTANT After the first processing, you might see an error message stating that the total number of results is too large for a single batch (more than 500 000). Return to the Processing pages in the method (refer to Chapter 6, “Using the Method Development Mode for Unknown Screening Methods,” in the *TraceFinder Lab Director User Guide*) and make one or more of these parameter adjustments:

- Limit the RT range.
- Shorten the signal range.
- Specify a lower value for the Number of Top Matches.
- Specify a Simple Search instead of an Exhaustive Search.
- Specify Top Peaks instead of All Peaks.
- Limit the number of search types.

Figure 117. Submit Options dialog box

Table 65. Submit Options dialog box parameters (Sheet 1 of 3)

Parameter	Description
User Name	Name of the current user.
Samples	Number of samples to be submitted for acquisition, processing, or reporting.
Workflow Steps to Perform	
Acquire Data	<p>Submits the current batch to acquisition.</p> <ul style="list-style-type: none"> When all submitted samples have been previously acquired, this option is (by default) not selected. When one or more samples in the batch have not been acquired, this option is (by default) selected.

Table 65. Submit Options dialog box parameters (Sheet 2 of 3)

Parameter	Description
Process Data	<p>Processes the data for the current batch using any of the following options:</p> <p>Peak Detect: Performs peak detection for all method types. You can process the data with or without performing peak detection. For example, you might want to turn off peak detection when reprocessing samples.</p> <p>Identify: Performs identification for unknown screening methods.</p> <p>With RT Alignment: Performs retention time alignment for unknown screening methods. This produces the heat map and group averages data in the Unknown Screening View. When you select this option, the application automatically selects the Peak Detect option.</p>
	<p>Note You can use Unknown Screening features in conjunction with Quantitation or Target Screening methods.</p> <p>For specifying processing data for batches that also use Quantitation methods, see To submit samples in the batch.</p> <p>For specifying processing data for batches that also use Target Screening methods, see To submit samples in the batch.</p>
Create Reports	Creates reports for the current batch.
Submit as a Priority Sequence?	
Priority Sequence	<p>With multiplexing activated, places the batch immediately after the currently acquiring batch.</p> <p>Without multiplexing activated, specifies one of the following priority options to place the batch in the queue:</p> <p>Next Available Batch: Places the batch immediately after the currently acquiring batch.</p> <p>Next Available Sample: Places the batch immediately after the currently acquiring sample.</p>
	<p>Note When you select Full Sequence Submission in the Configuration console, these options are unavailable because the current batch and the current sample are, in effect, the same thing.</p>

Table 65. Submit Options dialog box parameters (Sheet 3 of 3)

Parameter	Description
Acquisition pane	
Device Name	Lists all configured instruments. If the instrument that you want to use is not configured, close the TraceFinder application, configure the instrument, and then reopen the application. You cannot configure an instrument while the TraceFinder application is running. Available only when you select the Acquire Data check box.
Use	Specifies the instruments used for this acquisition. Available only when you select the Acquire Data check box.
Start Device	Specifies the instrument that initiates the communication with the other instruments. This is usually the autosampler. Available only when you select the Acquire Data check box.
Start When Ready	Starts the specified device when all the instruments are ready to acquire data. When this is cleared, individual instruments can start at different times and then must wait for the last instrument to be ready.
Post-run System State	Specifies the system state after it acquires the last batch: On (default), Standby, or Off.
Buttons	
Hide/Show Details	Collapses or expands the acquisition details of the Submit Options dialog box.
OK	Begins the selected processes.
Cancel	Closes the Submit Options dialog box without submitting any tasks.

❖ **To view the output files**

Locate the files to view from the following directories:

The application writes saved batches to the project folder:

...\\TraceFinderData\\4.0\\Projects\\...

For each acquired sample, the application writes an RSX file to the batch Data folder:

...\\TraceFinderData\\4.0\\Projects\\...\\Data

The application saves method information to the batch Methods folder:

...\\TraceFinderData\\4.0\\Projects\\...\\Methods

The application writes the reports to the batch Reports folder:

...\\TraceFinderData\\4.0\\Projects\\...*batch*\\Reports

Saving a Batch to a New Location

You can move the current batch to a different project folder, or you can make a copy of the current batch and save the copy to a different project folder.

Follow these procedures:

- [To save a batch to another project folder](#)
- [To move a batch to another folder](#)
- [To create new project folder](#)
- [To delete project folders](#)
- [To rename project folders](#)

❖ To save a batch to another project folder

1. Choose **File > Save > Save Batch As** from the Batch View main menu in the Analysis mode.

The [Save Batch As Dialog Box](#) opens.

2. Select a storage location.

The default storage location is C:\TraceFinderData\4.0\Projects.

3. Select or create a project folder.

4. Type a name for the new batch.

If you are saving the batch to a different folder, you must give it a unique name. You cannot overwrite an existing batch in a folder.

5. Click **Save**.

When you save the batch to a different folder, the reports reflect the original project folders and the application does not save the calibration history.

❖ To move a batch to another folder

1. Choose **File > Save > Move Batch** from the Batch View main menu in the Analysis mode.

The [Save Batch As Dialog Box](#) opens.

2. Select a storage location.

The default storage location is C:\TraceFinderData\4.0\Projects.

3. Select or create a project folder.


4. Type a name for the new batch.

You must give the batch a unique name in the new subproject folder. You cannot overwrite an existing batch.

5. Click **Save**.

When you move the batch, the reports reflect the original project and subproject folders and the application does not save the calibration history.

❖ To create new project folder

1. In the Save Batch As dialog box, select the folder for which you will create a new lower-level folder.
 - You can select the main TraceFinderData\4.0\Projects folder and create a new folder under it.
 - You can select one of the existing folders and create a lower-level folder under it.
2. Click the **Create Folder** icon, .


The application adds a new lower-level folder to the selected folder.

3. Select the new folder name and type a name for the folder.

Folder names can contain spaces and special characters, except for the following special characters: \ / : + ? " ' < >

Note After you add a lower-level folder, you cannot rename the parent folder.


❖ To delete project folders

1. In the Save Batch As dialog box, select the folder to delete.
2. Click the **Delete Folder** icon, .

With no confirmation prompt, the application immediately removes the selected folder.

Note This feature is not available for folders that contain lower-level project or batch folders; you must first delete the lower-level project or batch folders.

❖ To rename project folders

1. In the Save Batch As dialog box, select the folder to rename.
2. Click the **Rename Folder** icon, .

Note This feature is not available for folders that contain lower-level project or batch folders; you must first delete the lower-level project or batch folders.

3. Type a new name for the folder and press ENTER.

The application saves the new folder name.

Save Batch As Dialog Box

Use the features in the Save Batch As dialog box to save a batch to a new name or to move a batch to a different project folder.

Figure 118. Save Batch As dialog box

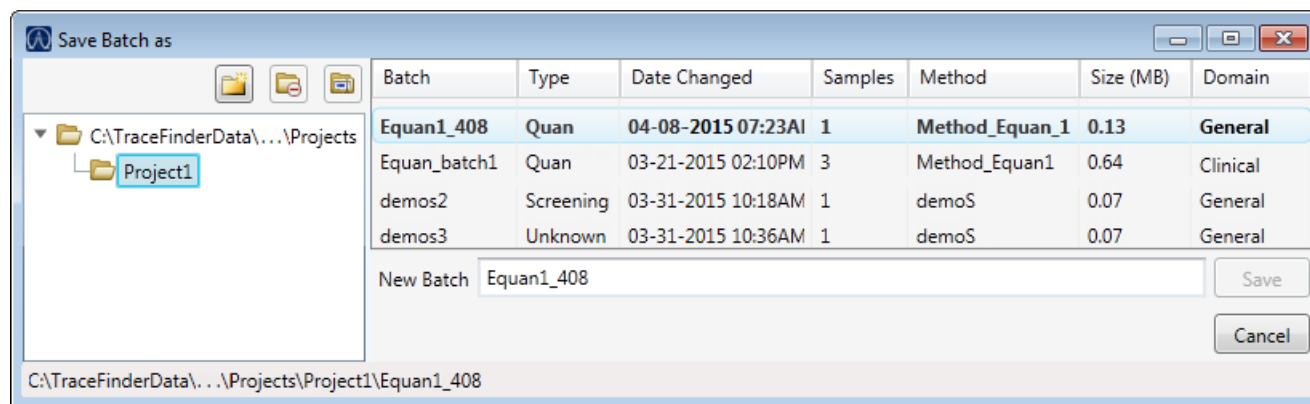


Table 66. Save Batch As dialog box parameters (Sheet 1 of 2)




Parameter	Description
 Create New Folder	<p>Adds one of the following:</p> <ul style="list-style-type: none"> When a drive is selected, adds a new project-level folder to the drive. When a project folder is selected, adds a subproject-level folder to the selected project. When a subproject folder is selected, adds a lower-level folder to the subproject. <p>Or, you can right-click and choose Create Folder.</p>
 Delete Folder	<p>With no confirmation prompt, immediately removes the selected folder.</p> <p>You cannot delete a folder that contains lower-level project or batch folders; you must first delete the lower-level project or batch folders.</p> <p>Or, you can right-click and choose Delete.</p>
 Rename Folder	<p>Renames the selected folder.</p> <p>You cannot rename a folder that contains lower-level project or batch folders; you must first delete the lower-level project or batch folders.</p> <p>Or, you can right-click and choose Rename.</p>
Batch table	
Batch	Name of batches in the selected project.
Type	Type of batch: Quan, Screening, or Unknown Only.
Date Changed	Date that the batch was last updated.
Samples	Number of samples in the batch.
Method	Name of the method used to create the batch.
Size	Size of the batch in megabytes.
Domain	TraceFinder domain in which the batch was created.

Table 66. Save Batch As dialog box parameters (Sheet 2 of 2)

Parameter	Description
New batch parameters	
New Batch	Name of the new batch to create. Note If the Create button is not activated, you have entered a name that is already used or you have not selected a method.
<i>Path</i>	Path to the project in the TraceFinderData\4.0\Projects folder where the batch is created.
Buttons	
Save	Saves the batch to the specified name and folder and opens the Batch View for the new batch.
Cancel	Closes the Save Batch As dialog box without saving the batch.
Shortcut menu commands	
Create Folder	Adds one of the following: <ul style="list-style-type: none"> • When a drive is selected, adds a new project-level folder to the drive. • When a project folder is selected, adds a subproject-level folder to the selected project. • When a subproject folder is selected, adds a lower-level folder to the subproject.
Delete Folder	Immediately removes the selected folder. There is no prompt to confirm that you want to delete the selected folder. You cannot delete a folder that contains lower-level project or batch folders; you must first delete the lower-level project or batch folders.
Rename Folder	Renames the selected folder. You cannot rename a folder that contains lower-level project or batch folders; you must first delete the lower-level project or batch folders.
Expand Child Nodes	Expands all project and subproject folders in the Project tree.
Collapse Child Nodes	Collapses all project and subproject folders in the Project tree.

Working in Data Review for Unknown Screening Batches

The Data Review view displays the data generated by the master method. Use Data Review to verify the data for identified peaks before you generate reports.

When you run samples in an experiment, you can check for peaks that are not already identified in either an associated target screening or quantitation method and take appropriate steps when the application finds these peaks and proposes whether they match one or more compounds using various identification sources.

IMPORTANT When unknown screening features are associated with a quantitation or target screening method, the unknown screening results do not include any compounds (other than internal standards) that were identified by the quantitation or target screening methods.

Follow these procedures:

- [To open the Data Review view](#)
- [To display or hide a pane on the Unknown Screening View page](#)
- [To move, dock, or float Data Review panes](#)
- [To restore the default layout](#)

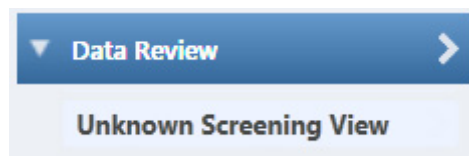
❖ To open the Data Review view

1. Click **Analysis** in the navigation pane.

The Analysis navigation pane opens.

2. Click **Data Review**.

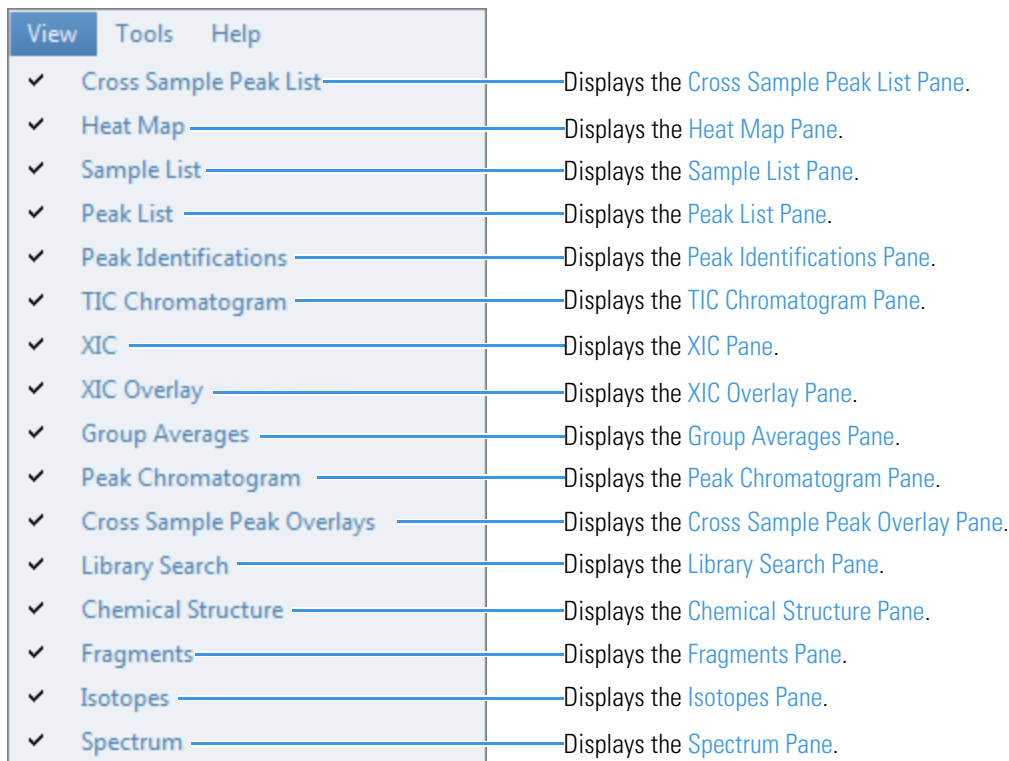
The Data Review navigation pane opens.



❖ **To display or hide a pane on the Unknown Screening View page**

From the View menu, choose to display or hide any of the following panes.

A check mark indicates a displayed pane.

❖ **To move, dock, or float Data Review panes**

Follow the instructions in [Appendix C, “Moving Data Review Panes.”](#)

❖ **To restore the default layout**

Choose **View > Restore Default Layout**.

The Unknown Screening View displays the default panes in their default locations.

Cross Sample Peak List Pane

Use the Cross Sample Peak List pane to compare peak values across all samples in a batch.

❖ **To use the Cross Sample Peak List pane**

1. In the Peak List pane, select a peak.

The application scrolls the Cross Sample Peak List pane to the specified peak and selects the peak.

2. Use the scroll bar at the bottom of the pane to scroll across the values for the selected peak in all samples and groups in the batch.

Figure 119. Cross Sample Peak List pane

Cross Sample Peak List						
	Selected	Retention Time	M/Z	Mass	Mono Isotopic Mass	Maximum Fold
	Aa ▼	= ▼	= ▼	= ▼	= ▼	= ▼
2	<input type="checkbox"/>	0.80	203.05	202.05	202.05	0.00
3	<input type="checkbox"/>	5.92	169.12	168.11	168.12	0.00
4	<input type="checkbox"/>	6.72	245.08	244.07	244.07	0.00

Apple_I	MS Area	Apple_I	Avg Area	Apple_I	CV	Apple_I	Fold
=	▼	=	▼	=	▼	=	▼
	0.00		1.00		0.00		0.00
	0.00		1.00		0.00		0.00
	157,846.41		157,846.41		0.00		0.00

Table 67. Cross Sample Peak List pane columns (Sheet 1 of 2)


Column	Description
Selected	Identifies peaks to display when you select the Selected option in the XIC Overlay pane, or identifies which peaks to process when you click the Submit ... for Processing icon,  . Do not confuse a check mark in the Selected column with the currently selected peak. The row for a currently selected peak is highlighted in blue.
Retention Time	The time after injection when the compound elutes. The total time that the compound is retained on the column.
M/Z	Mass-to-charge ratio found in the spectra for the peak. Assumes that the charge is 1.
Mass	The uncharged mass found in the spectra for the peak.
Monoisotopic Mass	The most abundant mass in the spectra for the peak.
Maximum Fold	The maximum difference of the group average response from the control average response.
Compound Name	Displays the compound name found in the library or compound database. This column is displayed only when the peaks are aligned and you have performed a simple search. Refer to Chapter 6, “Using the Method Development Mode for Unknown Screening Methods,” in the <i>TraceFinder Lab Director User Guide</i> .
Sample <i>n</i> MS Area	The MS peak area for each sample.
Default Average Area	The averaged response of all peak occurrences in the sample.

Table 67. Cross Sample Peak List pane columns (Sheet 2 of 2)

Column	Description
Default % CV	<p>The coefficient of variance for each sample.</p> <p>The coefficient of variance is calculated as the standard deviation of the multiple samples of one level, multiplied by 100, and divided by the average of the multiple samples of that level. This calculation is based on the areas of the peaks.</p>
Default Fold	The maximum difference of the sample average response from the control average response.
<i>Group n</i> Average Area	The averaged response of all peak occurrences in the group. The minimum value is 1, even when the peak never occurs in the group.
<i>Group n</i> % CV	<p>The coefficient of variance for each sample.</p> <p>The coefficient of variance is calculated as the standard deviation of the multiple samples of one level, multiplied by 100, and divided by the average of the multiple samples of that level. This calculation is based on the areas of all the peaks in the group of samples.</p>
<i>Group n</i> Fold	The maximum difference of the average response of all samples in a group from the control average response.

Heat Map Pane

Use the Heat Map pane to display the response of each peak occurrence in the batch. The Heat Map pane displays all MS Area values for all peaks in all samples in the batch that are above the peak threshold value specified in the method. When you select a peak in the Heat Map pane, the application displays the associated results for the selected peak in all panes of the Unknown Screening view.

The Heat Map pane is available only when you select the Alignment and Gap Filling option in the method (refer to Chapter 6, “Using the Method Development Mode for Unknown Screening Methods,” in the *TraceFinder Lab Director User Guide*) or select the RT Alignment option when you submit the batch (see [Submitting a Batch](#)).

Follow these procedures:

- [To use the Heat Map pane](#)
- [To use the Heat Map pane with the TIC Chromatogram pane](#)
- [To use the Heat Map pane with the XIC chromatogram pane](#)

❖ To use the Heat Map pane

1. Click the **Heat Map** tab to display the pane.

For detailed descriptions of all parameters in the Heat Map pane, see [Heat Map pane columns](#).

2. In the Peak List pane, select a peak.

The application scrolls the Heat Map pane to the specified peak and selects the peak.

For each found peak in the batch, the application assigns a color code indicating the relative MS area value from highest (orange) to lowest (green).

Highest	5,054,199.00
	2,041,157.25
	388,806.75
	220,048.25
	148,246.16
Lowest	102,254.38

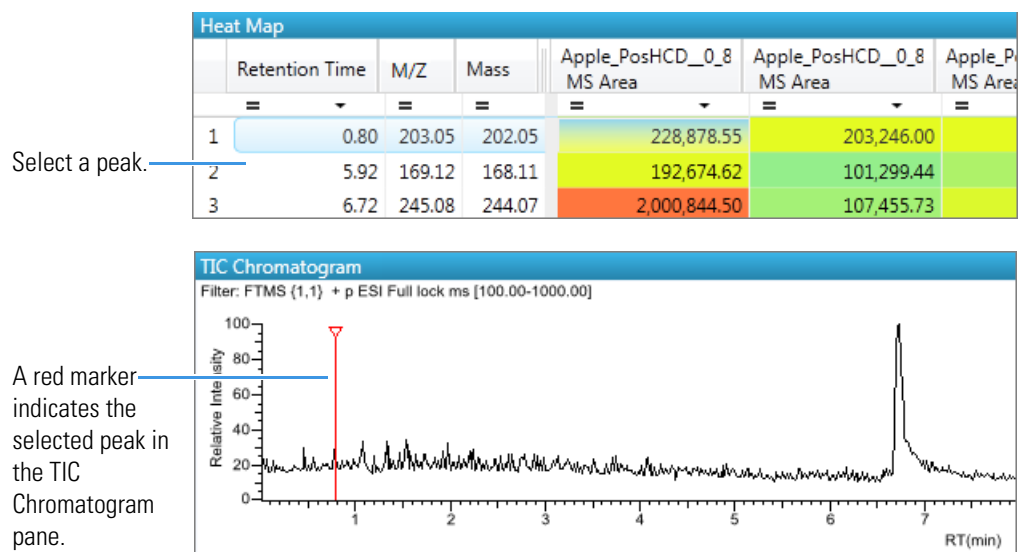
❖ To use the Heat Map pane with the TIC Chromatogram pane

1. Click the **TIC Chromatogram** tab to display the pane.

The total ion current (TIC) chromatogram pane displays the relative intensity of all found peaks in the batch.

2. In the Heat Map pane, select a peak.

The application identifies the selected peak in the TIC Chromatogram pane and indicates it with a red marker.



Or, you can select a peak in the TIC Chromatogram pane, and the application scrolls the heat map to the selected peak.

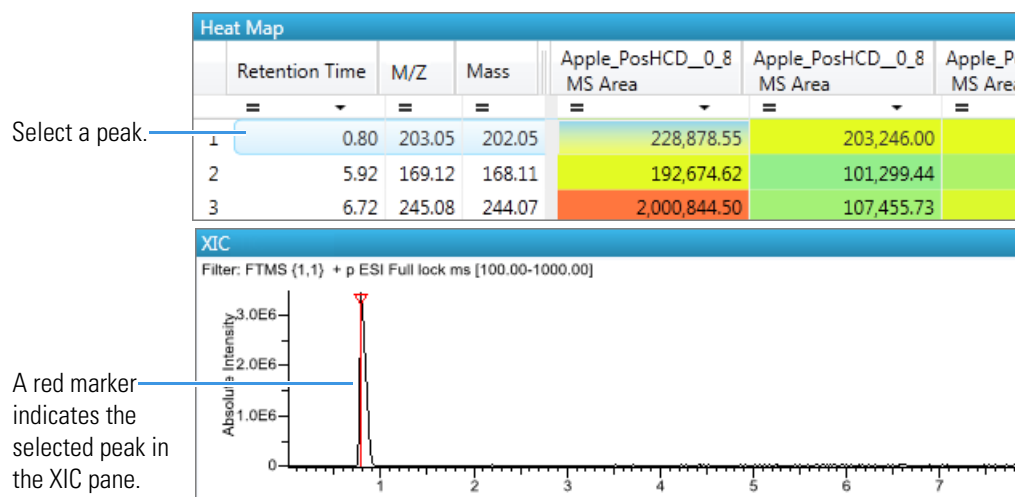
❖ **To use the Heat Map pane with the XIC chromatogram pane**

1. Click the **XIC** tab to display the pane.

The extracted ion chromatogram (XIC) pane displays the absolute intensity of the selected peak.

2. In the Heat Map pane, select a peak.

The application identifies the selected peak in the XIC, scales the chromatogram's *y* axis to fit the peak, and indicates it with a red marker.



Or, you can select a peak in the XIC pane, and the application scales the chromatogram to fit the peak, indicates it with a red marker, and scrolls the heat map to the selected peak.

Figure 120. Heat Map pane

Heat Map						
	Retention Time	M/Z	Mass	Apple_PosHCD_0_81_01 MS Area	Apple_PosHCD_0_8 MS Area	Apple_PosHCD_4_0 MS Area
1	5.92	169.12	168.11	192,674.62	101,299.44	112,746.64
2	6.72	245.08	244.07	2,000,844.50	107,455.73	170,410.11
3	6.72	163.04	162.03	1,156,472.00	205,919.31	107,573.62

Table 68. Heat Map pane columns

Column	Description
Retention Time	The time after injection when the compound elutes. The total time that the compound is retained on the column.
M/Z	Mass-to-charge ratio found in the spectra for the peak. Assumes that the charge is 1.
Mass	The uncharged mass found in the spectra for the peak.
Sample <i>n</i> MS Area	The MS peak area for each sample.

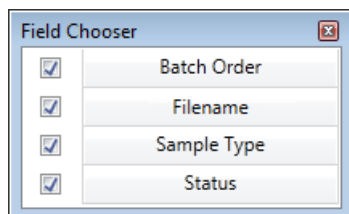
Sample List Pane

Use the [Sample List pane](#) to select a sample of interest. The application displays the associated results for the selected sample in all panes of the Unknown Screening view.

❖ To hide or display columns in the Sample List pane

1. Click the **Field Chooser** icon, , in the upper left corner of the pane.

The Field Chooser displays all available columns of data for the Sample List pane.



2. Select the check box for each column that you want to display, or clear the check box for each column that you want to hide.

The application immediately displays or hides the column in the Sample List pane.


3. When you are finished modifying the column display, click  to close the Field Chooser.

Figure 121. Sample List pane









Sample List				
	Batch Order	Status	Filename	Sample Type
	=	Aa	Aa	Aa
1	1		Unknown1	Specimen
2	2		Unknown2	Specimen
3	3		Unknown3	Specimen
4	4		Unknown4	Specimen

Table 69. Sample List pane columns (Sheet 1 of 2)

Column	Description
Batch Order	The order in which the samples were acquired.
Filename	A user-defined name that identifies a sample.

Table 69. Sample List pane columns (Sheet 2 of 2)

Column	Description
Sample Type	Defines how the application processes the sample data. Each sample is classified as one of the following sample types: Specimen, QC, Solvent, Calibrator, Hydrolysis, Unextracted, or Negative. Default: Specimen
Status	 Sample is not acquired.  Sample is acquired but not processed.  Sample is acquired and processed.  Sample is currently acquiring.

Peak List Pane

Use the [Peak List pane](#) to select a specific peak found in the selected sample. The Peak List displays each peak that the application identified in the sample.

❖ To display information for a peak

In the Peak List, click anywhere in the peak row.

The application displays information for the selected peak in the other Data Review panes.

Note Do not confuse a check mark in the Selected column with the currently selected peak. The row for the current, selected peak is highlighted in blue. A selected check box in the Selected column does not indicate the current, selected peak (see [Selected](#)).

Figure 122. Peak List pane

Peak List							
Selected	Peak ID	M/Z	Retention Time	Area	Height	Potential ID	Mass
1	peak @ 0.80 203.05	203.05	0.80	203,246.00	3,412,897.50	7	202.04
2	peak @ 11.48 312.3	312.36	11.48	455,449.00	7,354,474.69	7	311.54
3	peak @ 13.04 338.3	338.34	13.04	262,451.75	3,939,006.44	7	337.53
4	peak @ 14.20 278.9	278.97	14.20	166,002.72	513,232.62	9	277.96

Mono Isotopic Mass	Charge State	Filter String	NIST	mzVault
284.27	1	FTMS + p ESI Full ms [100.00-1000.00]	N/A	N/A
100.00	1	FTMS + p ESI Full ms [100.00-1000.00]	N/A	N/A
293.15	1	FTMS + p ESI Full ms [100.00-1000.00]	N/A	N/A
260.07	1	FTMS + p ESI Full ms [100.00-1000.00]	N/A	N/A

Elemental Composition	Database	Chemspider
C18H37O2	N/A	N/A
N/A	N/A	N/A
C13H23O4N2Na	N/A	N/A
C7H21O2N2S3	N/A	N/A

Table 70. Peak List pane columns (Sheet 1 of 2)


Column	Description
Selected	Identifies peaks to display when you select the Selected option in the XIC Overlay pane, or identifies which peaks to process when you click the Submit ... for Processing icon,  . Do not confuse a check mark in the Selected column with the currently selected peak. The row for the current, selected peak is highlighted in blue.
Peak ID	Identifies the peak as "peak @ <i>retention time m/z</i> ." Example: peak @ 0.80 203.05
<i>m/z</i>	Mass-to-charge ratio found in the spectra for the peak. Assumes the charge is 1.
Retention Time	The time after injection when the compound elutes. The total time (in minutes) that the compound is retained on the column.
Area	The area obtained by integrating peak intensities from the start to the end of the peak.
Height	The distance from the peak maximum to the peak base, measured perpendicular to the ordinate.
Potential ID	The number of returned peak IDs.
Mass	The uncharged mass found in the spectra for the peak.

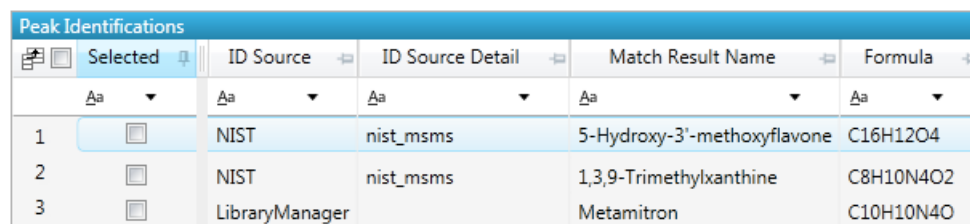
Table 70. Peak List pane columns (Sheet 2 of 2)

Column	Description
Monoisotopic Mass	The sum of the masses of the most abundant isotope.
Charge State	The charge state of the ion (the z value in m/z). For example, a charge state of 2 with a negative polarity means that the compound has 2 more electrons than protons. Valid range: 1 through 10 Default: 1
Filter String	Filter used to identify the peak. Specified in the raw data file or the master method.
NIST	The NIST match result name for the identified peak. If there is no match or if the Library search was not specified in the method, this shows as "N/A."
mzVault	The library match result name for the identified peak. If there is no match or if the Library Search was not specified in the method, this shows as "N/A."
Elemental Composition	The chemical formula for the identified peak. If there is no match or the Elemental Composition search was not specified in the method, this shows as "N/A."
Database	The compound database match result name for the identified peak. If there is no match or the Database Search was not specified in the method, this shows as "N/A."
ChemSpider	The ChemSpider database match result name for the identified peak. If there is no match or the ChemSpider Search was not specified in the method, this shows as "N/A."

Peak Identifications Pane

Use the Peak Identifications pane to display the name and formula for identified peaks and the source of the identification.

Figure 123. Peak Identifications pane



	Selected	ID Source	ID Source Detail	Match Result Name	Formula
1	<input checked="" type="checkbox"/>	NIST	nist_msms	5-Hydroxy-3'-methoxyflavone	C16H12O4
2	<input type="checkbox"/>	NIST	nist_msms	1,3,9-Trimethylxanthine	C8H10N4O2
3	<input type="checkbox"/>	LibraryManager		Metamitron	C10H10N4O

Table 71. Peak Identifications pane columns (Sheet 1 of 2)


Column	Description
Selected	Identifies peaks to display when you select the Selected option in the XIC Overlay pane, or identifies which peaks to process when you click the Submit ... for Processing icon,  . Do not confuse a check mark in the Selected column with the currently selected peak. The row for a currently selected peak is highlighted in blue.
ID Source	Specifies the search type used to identify the peak (Database Search, Library Search [NIST or mzVault], Elemental Composition, or ChemSpider Search). Refer to the instructions “To specify the search options” in Chapter 6, “Using the Method Development Mode for Unknown Screening Methods,” in the <i>TraceFinder Lab Director User Guide</i> .
ID Source Detail	Specifies the name of the library, the compound database, or the ChemSpider hyperlink where the peak was identified.
Match Result Name	Specifies the compound name that was found in the library, in the compound database, or in the ChemSpider hyperlink.
Formula	Specifies the chemical formula that was found in the library, in the compound database, in the elemental composition, or in the ChemSpider hyperlink.
Isotope Pattern Score %	Specifies the percentage of matched isotopes in the total number of isotopes.
Number of Isotopes Matched	Specifies the number of isotopes matched in the calculated isotope spectra relative to the total number of isotopes used in the score calculation, in the format “x of y”, where <ul style="list-style-type: none"> x = the number of isotopes matching the elemental composition used for the Isotope Pattern Score calculation. y = the total number of isotopes considered in the Isotope Pattern Score calculation. This is the number of isotope peaks expected to be above the spectral noise.

Table 71. Peak Identifications pane columns (Sheet 2 of 2)

Column	Description
Score	Specifies the score from the library fit. When the application finds a match in the library, in the compound database, or through the ChemSpider hyperlink, this column displays the highest score associated with the Match Result Name parameter. Valid range: 1 through 100%
SI/Dot Product	Specifies the search index method used to search the NIST library.
RSI/Rev Product	(Reverse search index) Specifies the method used to search the NIST library. A reverse search compares a library entry to an unknown compound (whereas a forward search compares the mass spectrum of an unknown compound to a mass spectral library entry).
Library Probability Percent	Specifies the probability ranking from the NIST library. When you use a reverse search, this value is "N/A."
Fragments	Specifies the number of fragments matched in the calculated spectra relative to the total number of expected fragments.

TIC Chromatogram Pane

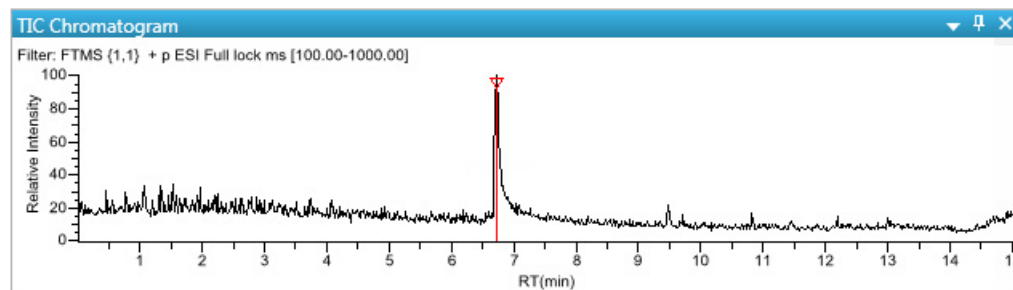
Use the TIC Chromatogram pane to display the relative intensity of a trace along the length of the sample data retention window.

❖ To focus on a specific peak

In the TIC Chromatogram pane, click any displayed peak.

The application indicates the selected peak with a red, vertical marker and updates the focus in the other Data Review panes with data for that peak.

Figure 124. TIC Chromatogram pane



XIC Pane

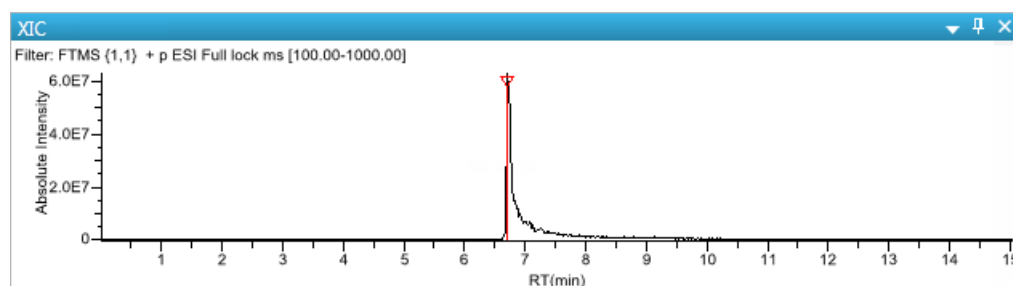
Use the XIC pane to display the absolute intensity of an extracted trace along the length of the sample data retention window.

❖ To focus on a specific peak

In the XIC pane, click any displayed peak.

The application indicates the selected peak with a red, vertical marker and updates the focus in the other Data Review panes with data for that peak.

Figure 125. XIC pane



XIC Overlay Pane

Use the XIC Overlay pane to view specific groups of peaks. You can choose to view all peaks, selected peaks only, or the top 20 most intense peaks (by area). The XIC Overlay plot is a collection of overlaid, extracted m/z ion plots that use a different color for each peak.

Follow these procedures:

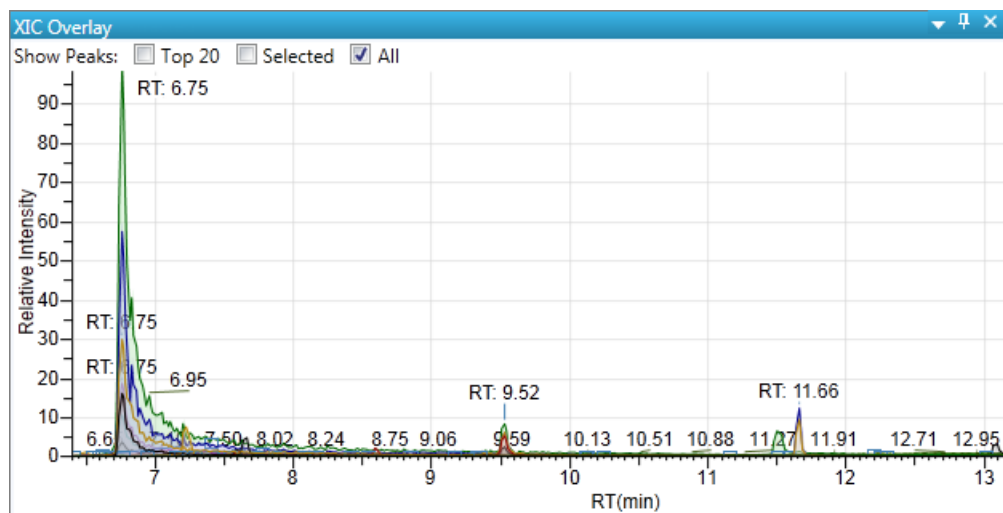
- To display all peaks in a sample
- To display only selected peaks in a sample
- To display the top 20 peaks in a sample
- To focus on a specific peak

❖ To display all peaks in a sample

1. In the Sample List pane, select the sample whose peaks you want to view.
2. In the XIC Overlay pane, select the **All** check box.

The application displays all the peaks in the selected sample. See [XIC Overlay plot showing all peaks](#).

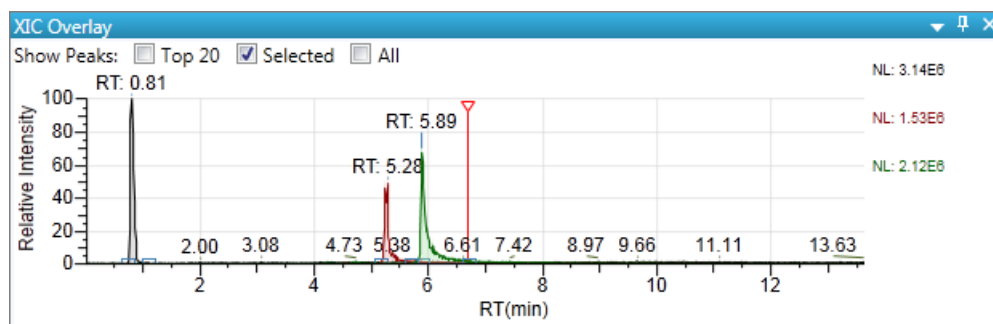
Note If there are more than 1000 peaks in the sample, performance might be slow.

Figure 126. XIC Overlay plot showing all peaks

❖ **To display only selected peaks in a sample**

1. In the Sample List pane, select the sample whose peaks you want to view.
2. In the Peak List pane, select the **Selected** check box for each peak that you want to view.
3. In the XIC Overlay pane, select the **Selected** option.

The application displays all selected peaks.

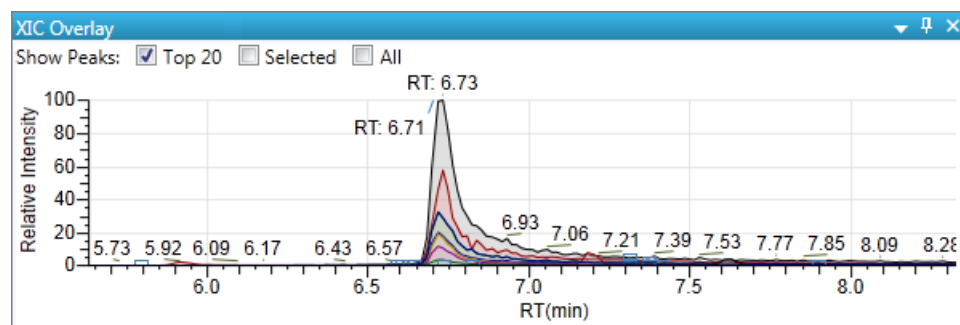
Figure 127. XIC Overlay plot showing selected peaks

❖ **To display the top 20 peaks in a sample**

1. In the Sample List pane, select the sample whose peaks you want to view.
2. In the XIC Overlay pane, select the **Top 20** check box.

The application displays the top 20 peaks in the selected sample.

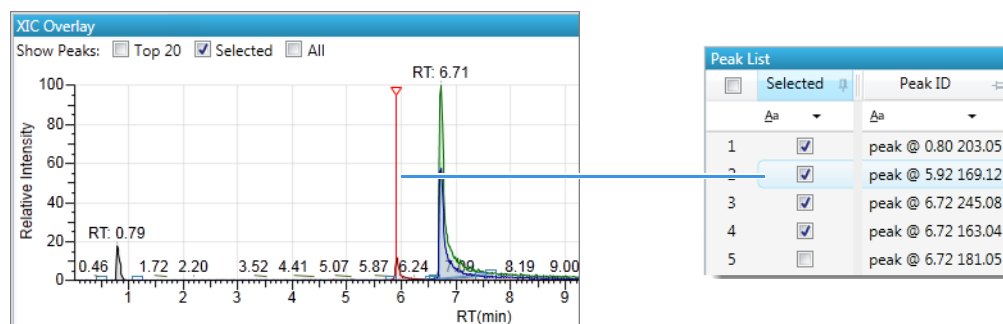
Figure 128. XIC Overlay plot showing the top 20 peaks



❖ To focus on a specific peak

In the XIC Overlay pane, click any displayed peak.

The application indicates the selected peak with a red, vertical marker and updates the focus in the other Data Review panes with data for that peak.



Group Averages Pane

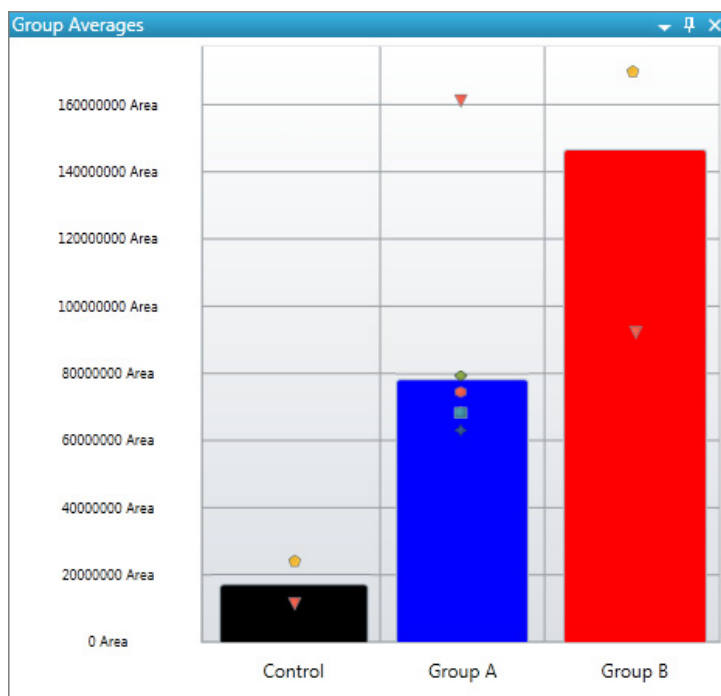
Use the Group Averages pane to compare the peak areas of different samples to a control group of samples.

The Group Averages pane is available only when you select the Alignment and Gap Filling option in the method (refer to Chapter 6, “Using the Method Development Mode for Unknown Screening Methods,” in the *TraceFinder Lab Director User Guide*) or select the RT Alignment option when you submit the batch (see [Submitting a Batch](#)).

The control group is displayed to the left, and other samples and groups are listed in the order in which they appear in the Batch View. To specify a group as the control group, see [Groups](#).

Each group displays a colored bar that indicates the average response of the group, as displayed in the Average Area column in the [Cross Sample Peak List Pane](#). The color of the bar matches the colors assigned to the samples and peaks in the [Cross Sample Peak Overlay Pane](#).

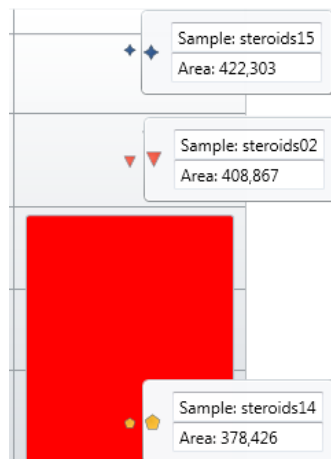
Figure 129. Group Averages pane



❖ **To display areas for each sample in the batch**

Hold the cursor over an indicator icon (▼, ◆).

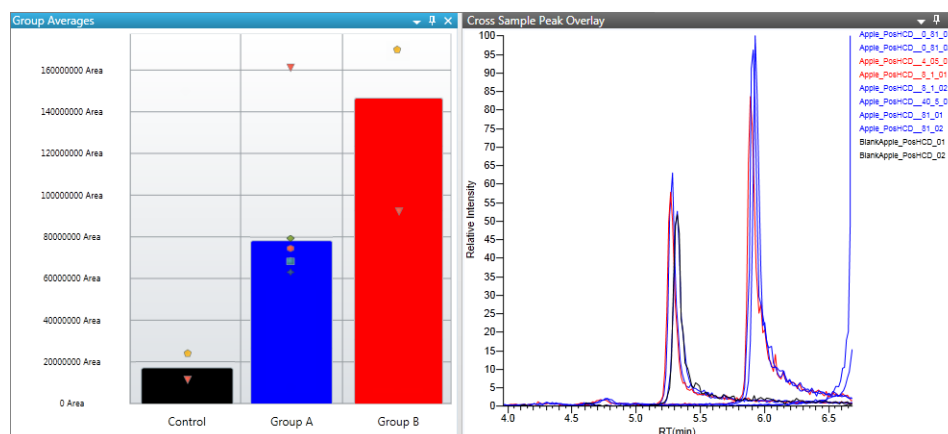
The application displays the area for the selected compound in each represented sample.



❖ To use the Group Averages pane with the Cross Sample Peak Overlay pane

Display both the Group Averages pane and the Cross Sample Peak Overlay pane.

The color of the bars in the Group Averages pane matches the colors assigned to the group samples and peaks in the Cross Sample Peak Overlay pane.



Peak Chromatogram Pane

Use the [Peak Chromatogram Pane](#) to display the selected chromatogram peak. The pane initially displays only the apex scan for the detected peak. You can manually integrate the peak and use the updated peak to generate reports.

Follow these procedures:

- [To zoom in on a peak](#)
- [To remove a peak](#)
- [To customize the labels in the display](#)
- [To display surrounding apex times](#)
- [To manually integrate a peak](#)
- [To switch between Method and Manual integration modes](#)

❖ To zoom in on a peak

1. In the chromatogram plot, drag the cursor to delineate a rectangle around the peak.

The delineated area expands to fill the view.

2. To restore the default view, right-click the chromatogram plot and choose **Reset Scaling**.

❖ To remove a peak

Right-click the Peak Chromatogram pane, and choose **Delete Peak**.

The application removes the peak displayed in the Peak Chromatogram pane. All data for this peak are removed from the Unknown Screening View panes.

❖ To customize the labels in the display

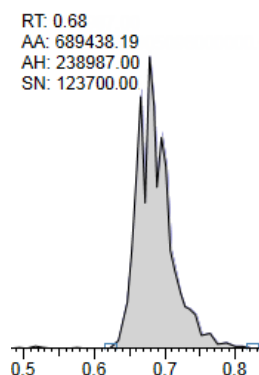
1. Right-click the Peak Chromatogram pane, and choose **Peak Display Settings**.

The [Peak Settings Dialog Box](#) opens.

2. To display labels for the peak retention time (RT), mass height (MH), mass area (MA), or signal-to-noise (SN), select the appropriate check box.
3. Click **OK** in the Peak Settings dialog box.

The chromatogram plot displays the selected labels.

Figure 130. Example of displayed labels

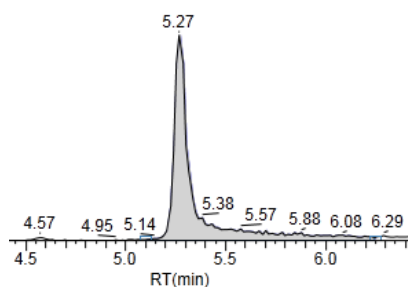
**❖ To display surrounding apex times**

1. Right-click the Peak Chromatogram pane, and choose **Peak Display Settings**.


The [Peak Settings Dialog Box](#) opens.

2. Select the **Show All Apex Times** check box.
3. Click **OK** in the Peak Settings dialog box.

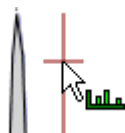
Figure 131. Example of displayed apex times



❖ To manually integrate a peak

1. Right-click the Peak Chromatogram pane and then choose **Swipe Manual Integration**, or click the **Swipe Manual Integration** icon, .

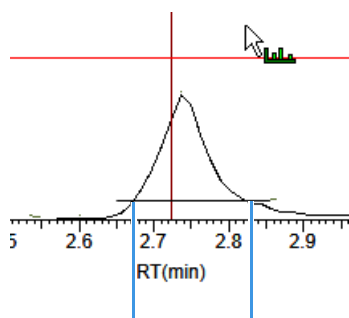
The cursor changes to indicate the manual integration mode.



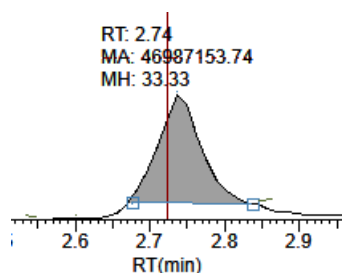
2. Drag the cursor to define the beginning and ending base points for the new peak. See [Manually integrated peak](#).

Note You must drag the cursor inside the x axis and y axis.

Figure 132. Manually integrated peak



Drag the cursor across the peak at the base points for the new peak.



The application identifies the peak and indicates the manual integration in the labels.

The application updates the area of the peak to the baseline that you just added. When you regenerate the reports, they display the manually modified data.

3. To return to the original peak as defined by the method, right-click and choose **Integration Settings > Method**.

❖ To switch between Method and Manual integration modes

Right-click the Peak Chromatogram pane and choose **Integration Settings > Manual** or **Integration Settings > Method**.

Initially, the method integration and manual integration settings that are stored for a compound and file are identical. When you switch modes, the saved result set does not change. However, when manual data are available, both the Peak Chromatogram plots and the result set update as you switch between method and manual modes.

As you switch between modes, each pane reflects the changes. The generated reports for the data identify the manual modifications.

Peak Chromatogram Pane

Use the Peak Chromatogram pane to display the selected chromatogram peak.

Figure 133. Peak Chromatogram pane

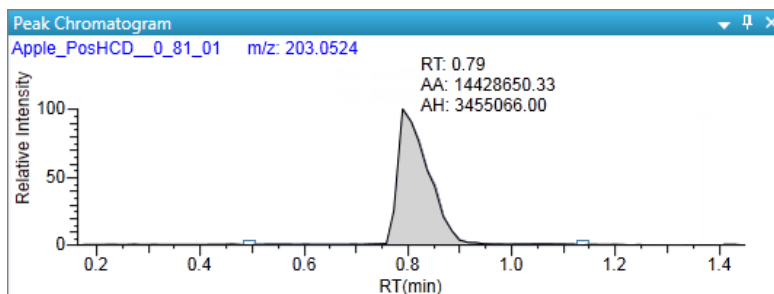




Table 72. Peak Chromatogram pane shortcut menu commands

Command	Description
Reset Scaling	Resets the original scaling after a zoom operation.
Copy to Clipboard	Copies the graphic display to the Clipboard.
Swipe Manual Integration	Switches to manual integration mode, where you can manually define the peak. Also available from the icon in the upper right corner,  .
Integration Settings (Manual or Method)	Switches the display between manual and method integration modes.
Delete Peak	Removes the peak and all data for this peak from the Unknown Screening View panes. Also available from the icon in the upper right corner,  .
Peak Display Settings	Opens the Peak Settings Dialog Box .

Peak Settings Dialog Box

Use the Peak Settings dialog box to custom the display settings.

Figure 134. Peak Settings dialog box

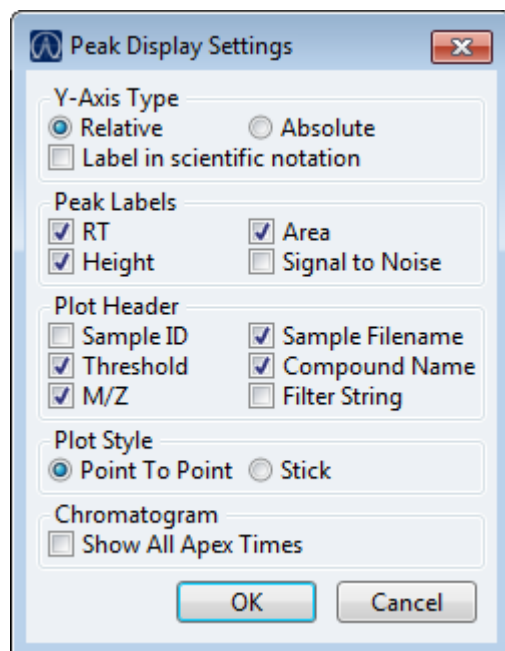
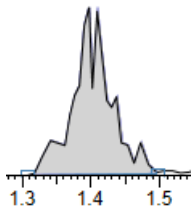
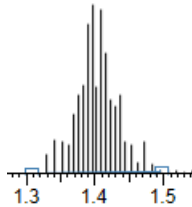


Table 73. Peak Settings dialog box parameters (Sheet 1 of 2)

Command	Description
Y-Axis Type	
Relative	Displays the y-axis scale from 0 through 100.
Absolute	Displays the y-axis scale from 0 to the actual value of the most intense peak in the group.
Label in Scientific Notation	Displays the y-axis scale in scientific notation.
Peak Labels	
RT	Displays the retention time (RT) label in the peak chromatogram pane.
Height	Displays the peak height (AH) label in the peak chromatogram pane.
Area	Displays the peak area (AA) label in the peak chromatogram pane.
Signal to Noise	Displays the signal-to-noise (SN) label in the peak chromatogram pane.
Plot Header	
Sample ID	Displays a user-defined, alphanumeric string that identifies a sample.
Threshold	Displays the peak threshold.

Table 73. Peak Settings dialog box parameters (Sheet 2 of 2)

Command	Description
<i>m/z</i>	Displays the mass-to-charge ratio found in the spectra for the peak. Assumes the charge is 1.
Sample Filename	Displays the name of the current, selected sample where the peak was found.
Compound Name	Displays the compound name match in the library or compound database.
Filter String	Displays the filter used to identify the peak. Specified in the raw data file or the master method.
Plot Style	
Point To Point	Displays chromatograms in the following format:
	
Stick	Displays chromatograms in the following format:
	
Chromatogram	
Show All Apex Times	Displays all surrounding apex times in the plot.

Cross Sample Peak Overlay Pane

Use the [Cross Sample Peak Overlay Pane](#) to compare instances of a selected peak across all samples in the batch. The application overlays all occurrences of the peak in the batch.

The application displays the names of all samples in the batch where the selected peak is found. Samples that are assigned to groups are color coded, and the peaks found in those samples are color coded in the plot. To assign samples to groups, see [Groups](#).

The retention time window displayed in the Cross Sample Peak Overlay pane is the RT window that you specified in the method.

Follow these procedures:

- To select the peak to display in the Cross Sample Peak Overlay pane
- To zoom in on a peak
- To use the Cross Sample Peak Overlay pane with the Group Averages pane

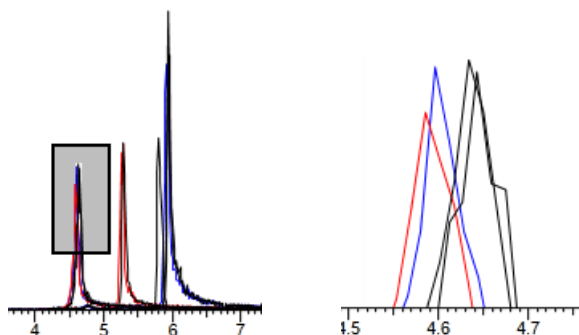
❖ To select the peak to display in the Cross Sample Peak Overlay pane

Select a peak in either the [Peak List Pane](#) or the [Cross Sample Peak List Pane](#).

❖ To zoom in on a peak

1. In the chromatogram plot, drag the cursor to delineate a rectangle around the peak.

The delineated area expands to fill the view.

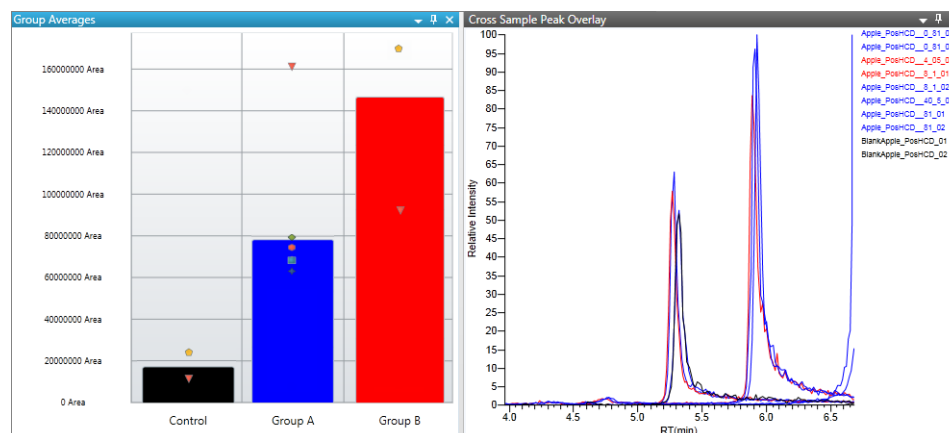


2. To restore the default view, right-click the chromatogram plot and choose **Reset Scaling**.

❖ To use the Cross Sample Peak Overlay pane with the Group Averages pane

Display both the Cross Sample Peak Overlay pane and the Group Averages pane.

The color of the bars in the Group Averages pane matches the colors assigned to the group samples and peaks in the Cross Sample Peak Overlay pane.



Cross Sample Peak Overlay Pane

Use the Cross Sample Peak Overlay pane to compare instances of a selected peak across all samples in the batch

Figure 135. Cross Sample Peak Overlay pane

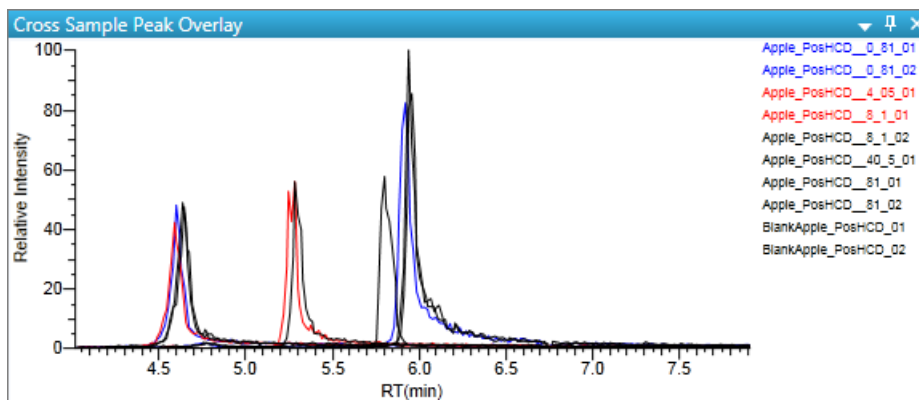


Table 74. Cross Sample Peak Overlay pane shortcut menu commands

Column	Description
Reset Scaling	Resets the original scaling after a zoom operation.
Copy to Clipboard	Copies the graphic display to the Clipboard.

Library Search Pane

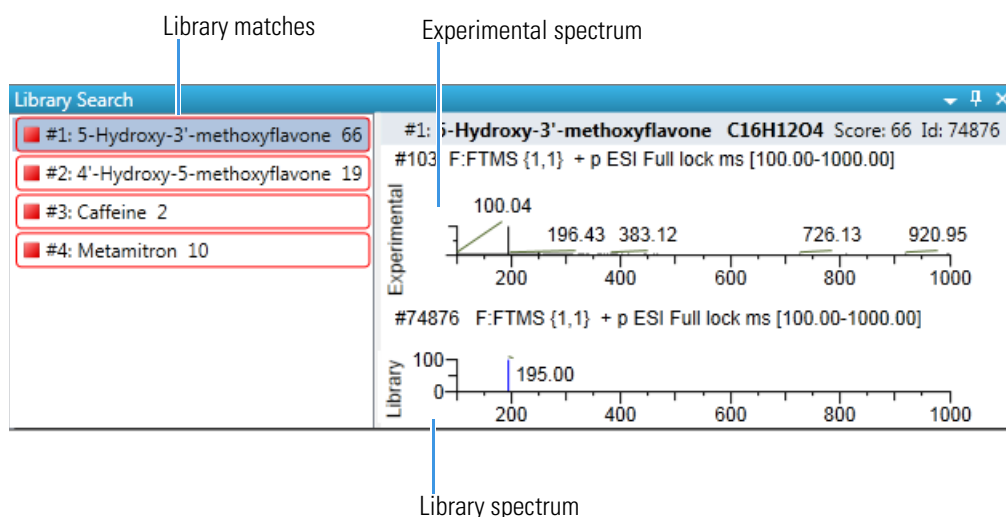
The [Library Search Pane](#) displays the best library matches for the selected peak, with the highest score listed first.

Follow these procedures:

- [To view the spectrum for a library match](#)
- [To display the difference between experimental and library spectra](#)

❖ To view the spectrum for a library match

In the Library Search pane, select a library entry.

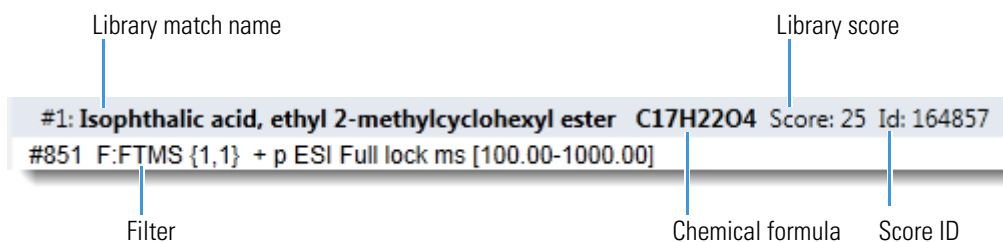


The Experimental plot displays the measured spectra for the selected compound. Library matches are colored red.

The Library plot displays the expected spectra in the library. Library matches are colored blue.

The spectra header displays the following:

- The library match name
- The chemical formula for the peak
- The score from the library fit when the library type is mzVault or NIST
- SI and RSI when the library type is NIST
- Dot Prod or Rev Dot Prod when the library is HRAM
- Molecular weight when the library type is NIST
- The ID record from the library when the library type is mzVault or NIST
- The scan number and scan filter when either is present in the library

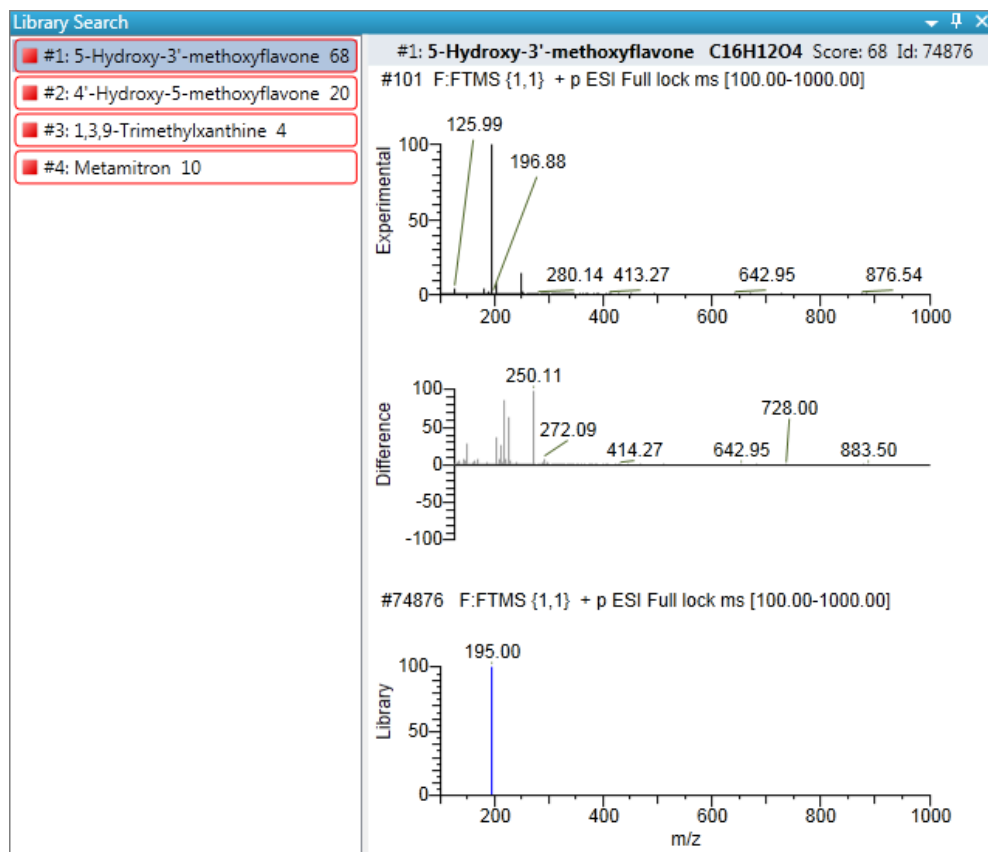


❖ To display the difference between experimental and library spectra

1. Right-click the spectrum plot and choose **Show Difference Plot**.

The application displays the Experimental and Library plots and a Difference plot that shows the difference between the experimental and library spectra. When you read a difference plot, the match between the experimental and library spectra is better when fewer peaks appear.

Figure 136. Library Search pane showing a Difference plot



2. To specify how the difference plot displays the mass tolerance, do the following:
 - a. Right-click the spectrum plot and choose **Set Difference Plot Mass Tolerance**.
 - b. In the dialog box that opens, select either **MMU** or **PPM** as the unit of measure.
The default value is the tolerance set in the NIST library or mzVault library.
 - c. Type a mass tolerance value.
 - d. Click **OK**.

Note This command is available only when you display the difference plot.

Library Search Pane

Use the Library Search pane to display the best library matches for the selected peak.

Figure 137. Library Search pane

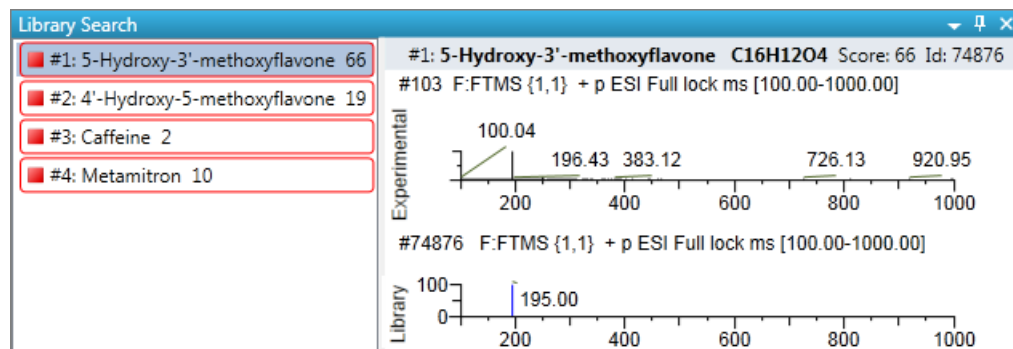


Table 75. Library Search pane shortcut menu commands

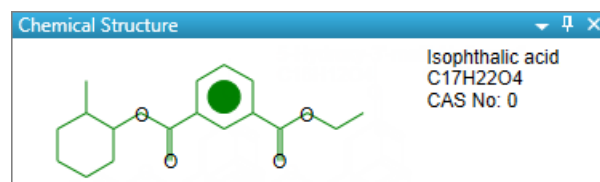
Command	Description
Reset Scaling	Resets the original scaling after a zoom operation.
Copy to Clipboard	Copies the graphic display to the Clipboard.
Display Overlay Spectra	Overlays the two spectrum displays, or stacks the Experimental and Library spectra.
Display Stack Spectra	When you display the Difference plot, this spectrum also displays the Difference spectrum.
Show/Hide Difference Plot	Displays or hides the Difference spectrum plot.
Set Difference Plot Mass Tolerance	Opens a dialog box where you specify the mass tolerance in either MMU or PPM units. Available only when you display the Difference Plot.

Chemical Structure Pane

Use the Chemical Structure pane to view the chemical formula and structure for the peak that is currently selected in the Peak List pane.

The application displays the chemical formula (and CAS number when available) that was found in the search database that you specified in the method. Refer to Chapter 6, “Using the Method Development Mode for Unknown Screening Methods,” in the *TraceFinder Lab Director User Guide*.

Figure 138. Chemical Structure pane



Fragments Pane

The Fragments pane displays the maximum number of fragments as specified in the unknown screening method. Refer to Chapter 6, “Using the Method Development Mode for Unknown Screening Methods,” in the *TraceFinder Lab Director User Guide*.

When a compound database search returns a match for a peak that has fragments, the Fragments pane displays the theoretical fragments.

When the data has an MS/MS element that is identified as belonging to the chromatographic peak, the Fragments pane displays the fragments found in the MS/MS scan.

The Fragments pane displays the fragments in one of two ways:

- [All Fragments](#)
- [Individual Fragments](#)

All Fragments

The All Fragments view displays a composite of all fragments found in the peak. The application displays the measured peak as a red line; the application displays the expected peak as a blue line. The application displays these headers for the All Fragments view:

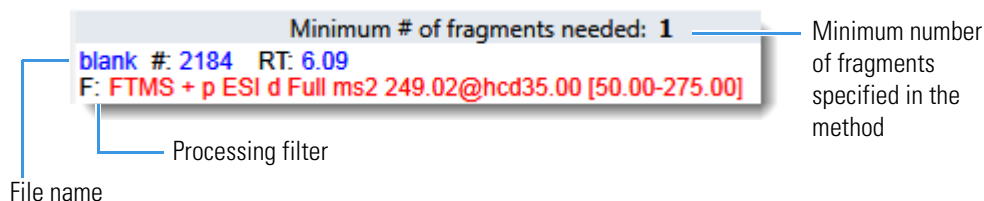


Figure 139. Fragments pane with overlaid spectra for all fragments

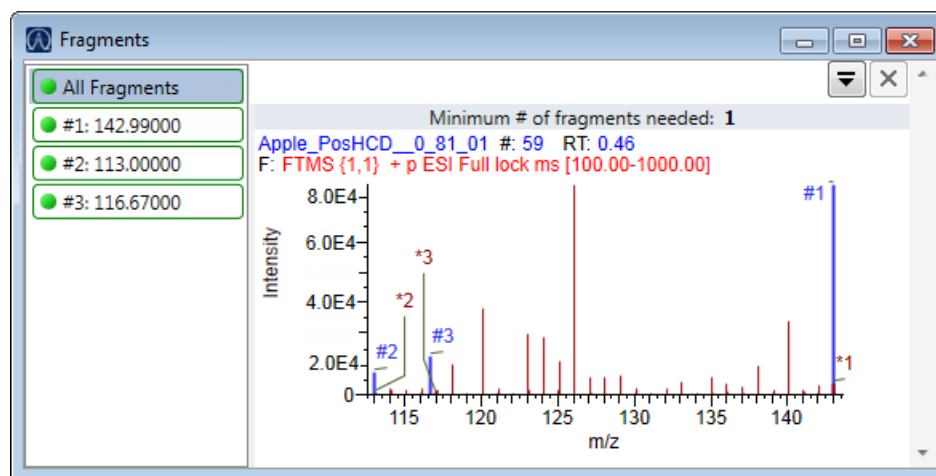
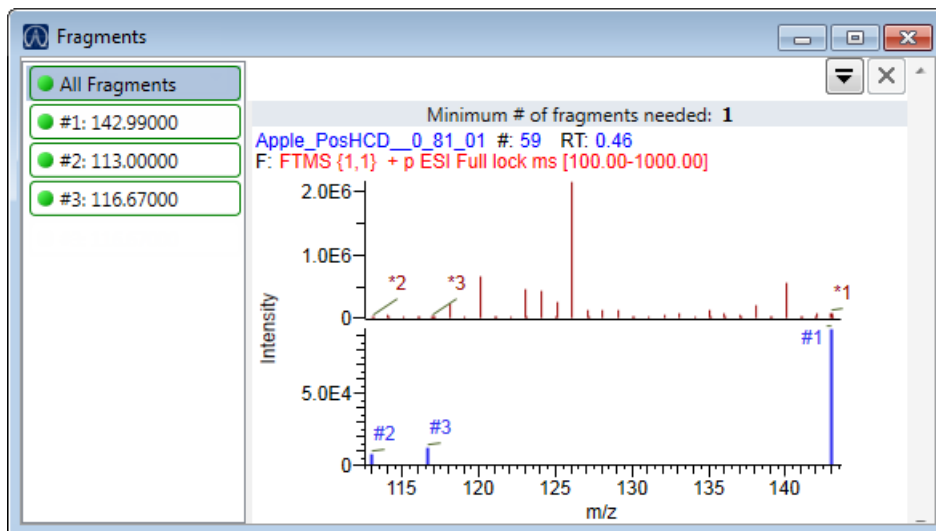


Figure 140. Fragments pane with stacked spectra for all fragments



Individual Fragments

The individual fragments view displays the expected and measured peaks for a single fragment. The application displays these headers for the individual fragments view:

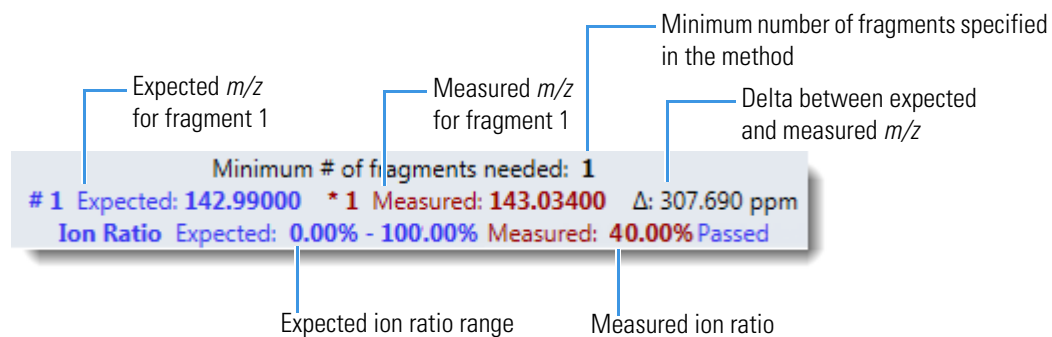
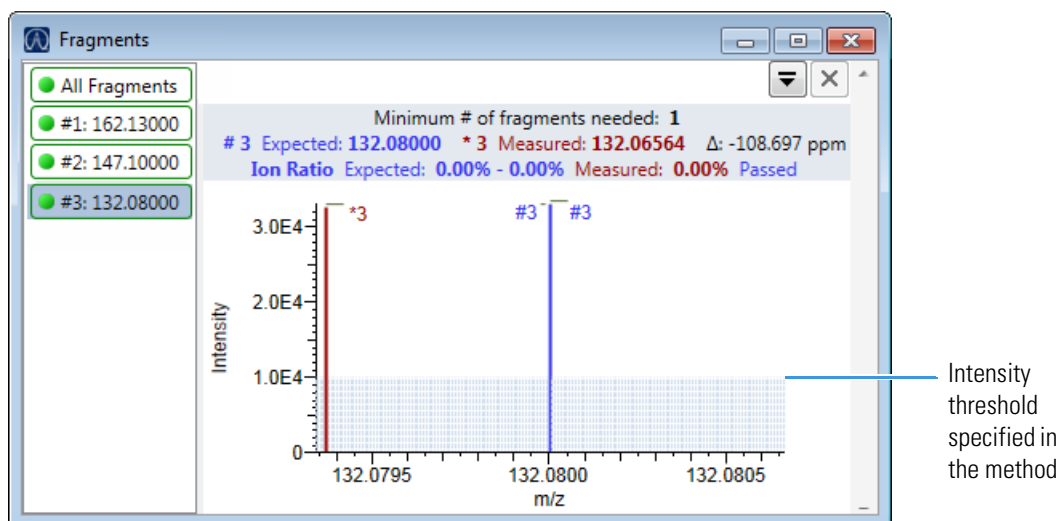
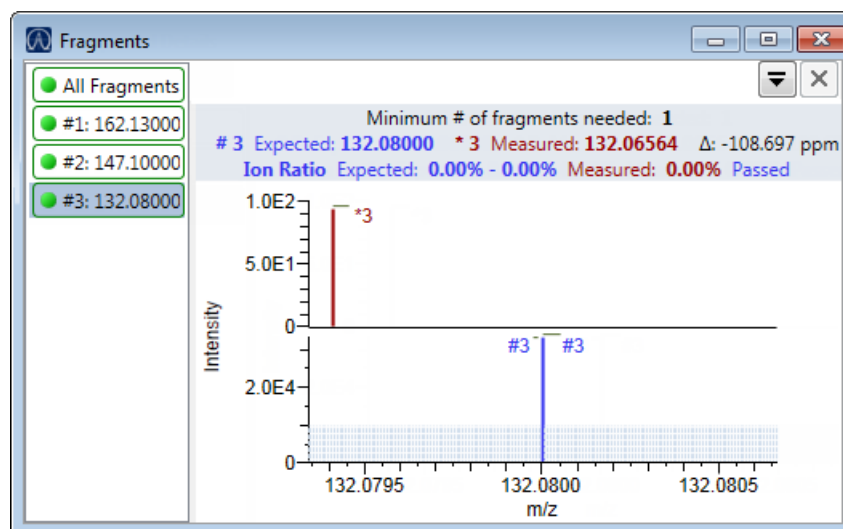


Figure 141. Fragments pane with overlaid spectra for a single fragment**Figure 142.** Fragments pane with stacked spectra for a single fragment

Isotopes Pane

The isotopes pane displays isotopic pattern results for all adducts of a compound according to the threshold and deviation parameters defined in the unknown method.

To identify or confirm the presence of a compound, the resulting score percentage from isotopic pattern matching must be higher than the specified fit threshold percentage.

- An isotope peak is not found if its intensity, relative to the monoisotopic ion's intensity, is more than the specified intensity deviation percentage away from the theoretical relative intensity of the isotope ion.
- An isotope peak is found if its measured m/z is less than the specified mass deviation amount away from its expected m/z .

To specify the peak threshold and deviation parameters, refer to Chapter 6, “Using the Method Development Mode for Unknown Screening Methods,” in the *TraceFinder Lab Director User Guide*.

The Isotopes pane displays the isotopes in one of three ways:

- [All Isotopes](#)
- [Multi-Isotopes](#)
- [Individual Isotopes](#)

All isotopes panes use a shortcut menu so that you can specify how to display the data. See [Isotopes pane shortcut menu commands](#).

All Isotopes

The All Isotopes view displays a composite of all isotopes found in the compound. The application scales the window with respect to the most intense isotope. The most intense isotope is usually the first isotope unless you are using halogenated compounds. The application displays the measured peak as a solid red line; the application displays the expected peak as a dashed blue line.

The application displays these headers for the All Isotopes view:

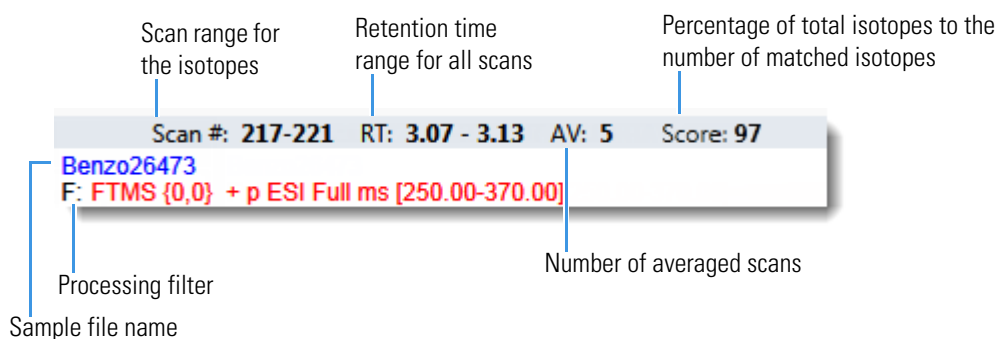


Figure 143. Isotopes pane with stacked spectra for all isotopes

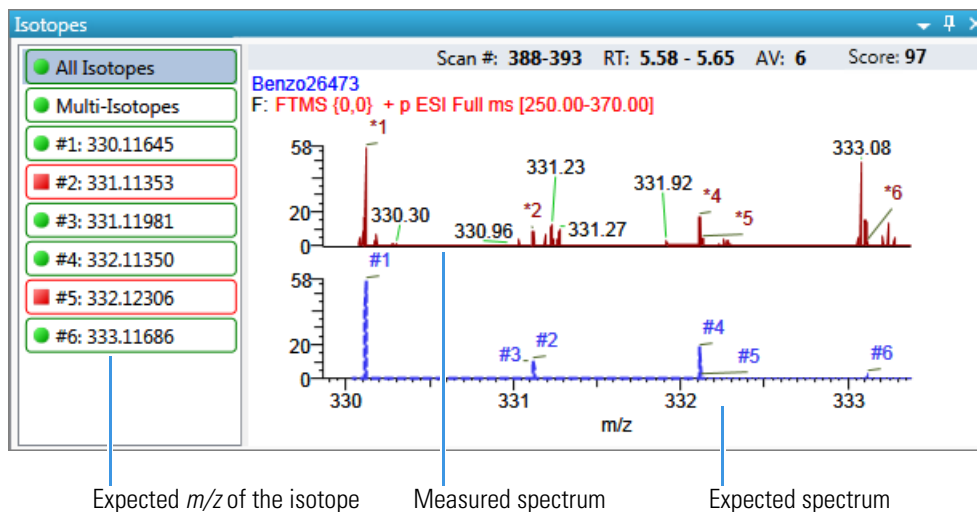
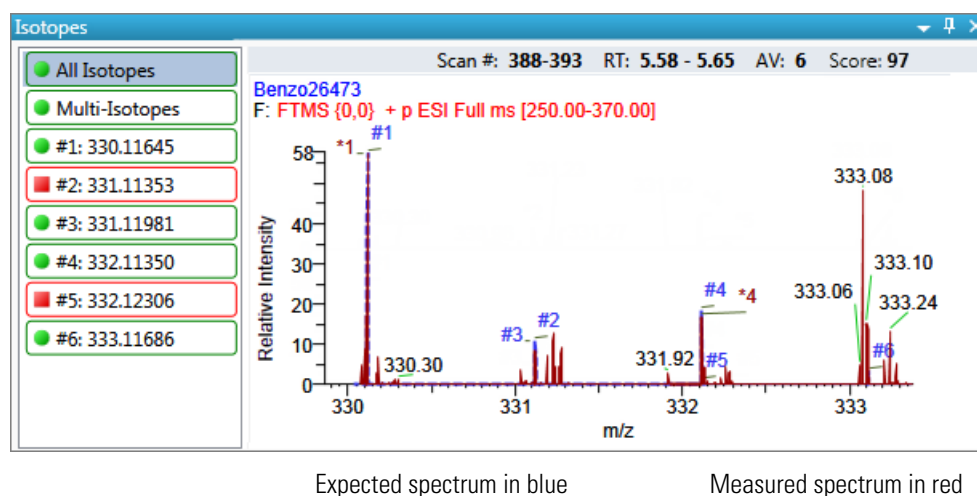


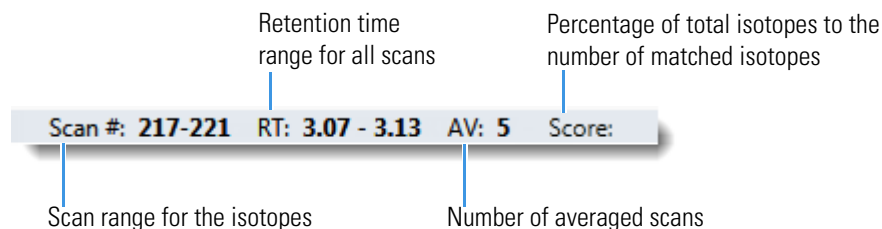
Figure 144. Isotopes pane with overlaid spectra for all isotopes



Multi-Isotopes

The Multi-Isotopes view displays individual plots for each isotope. You can individually stack or overlay the plots for each isotope.

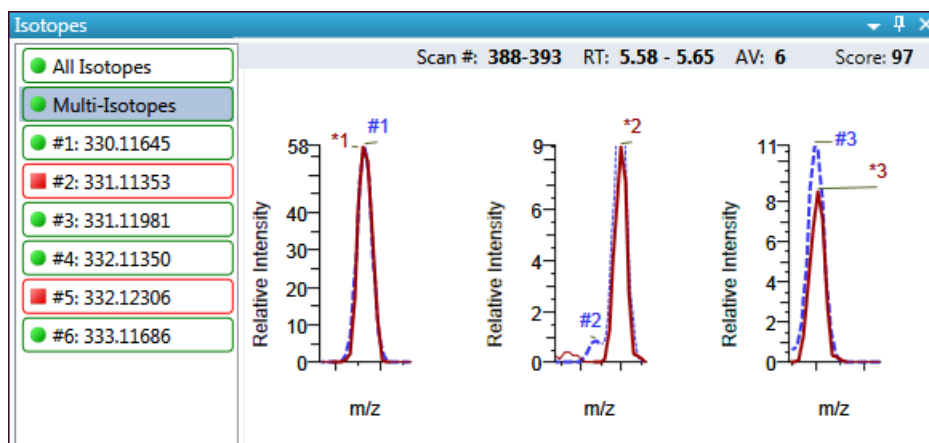
The application displays these headers for the Multi-Isotopes view:



6 Using the Analysis Mode for Unknown Screening Batches

Working in Data Review for Unknown Screening Batches

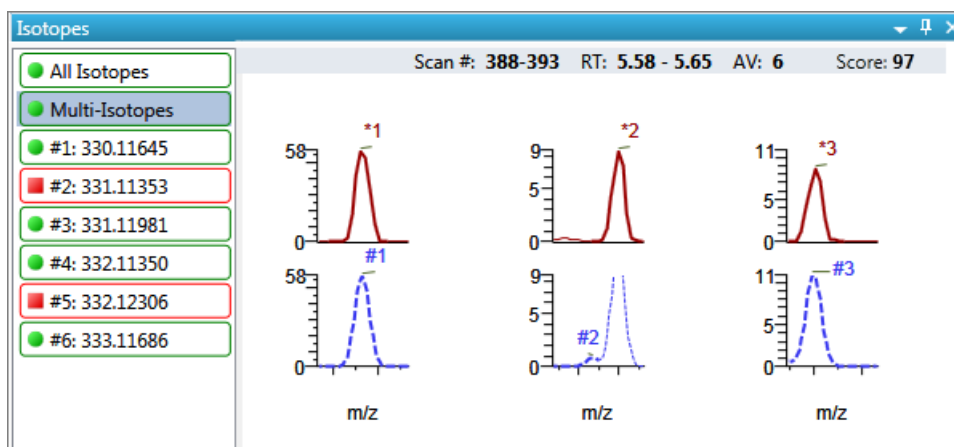
Figure 145. Isotopes pane with overlaid spectra for multi-isotopes



Expected spectrum in blue

Measured spectrum in red

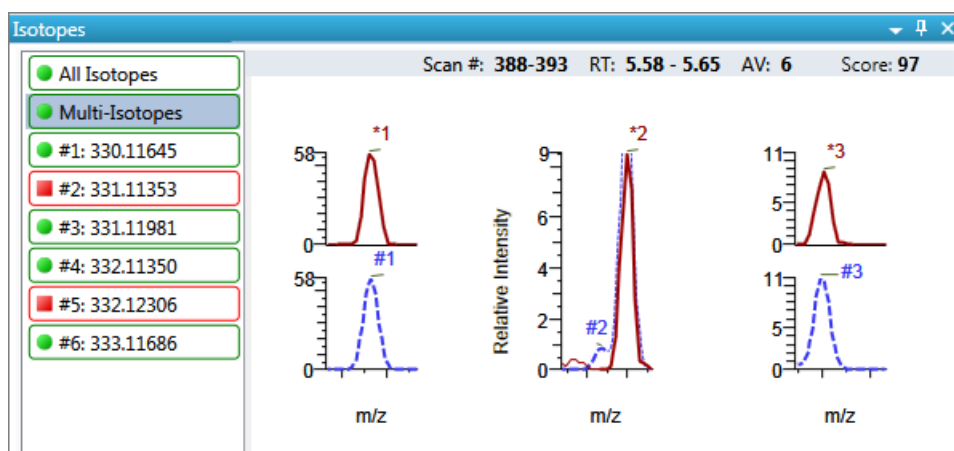
Figure 146. Isotopes pane with stacked spectra for multi-isotopes



Expected spectrum in blue

Measured spectrum in red

Figure 147. Isotopes pane with stacked and overlaid spectra for multi-isotopes



Expected spectrum in blue

Measured spectrum in red

Individual Isotopes

The individual isotopes view displays the expected and measured peaks for a single isotope.

The application displays these headers for the individual isotopes view:

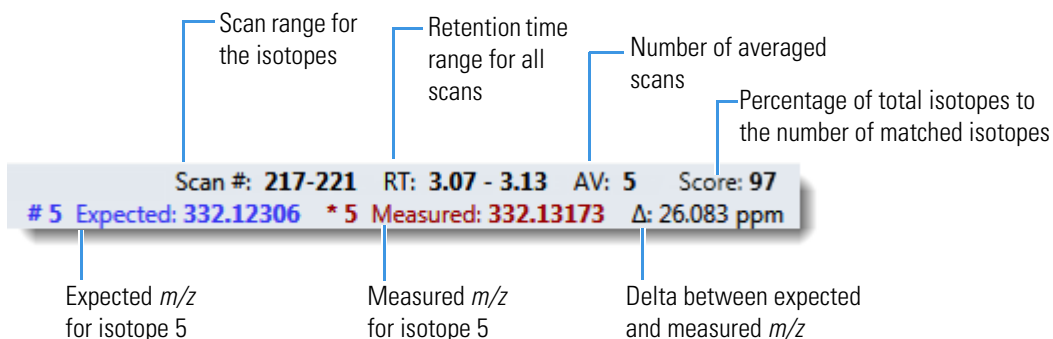


Figure 148. Isotopes pane with overlaid spectra for a single isotope

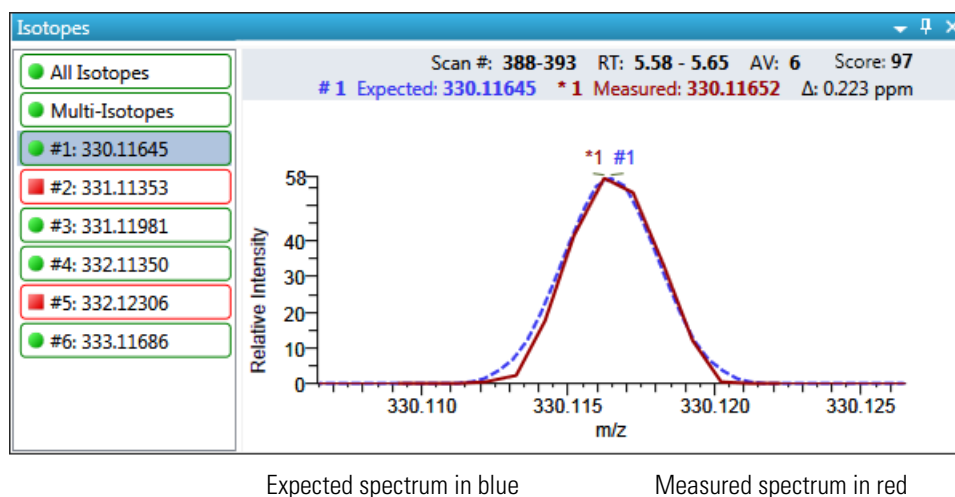


Figure 149. Isotopes pane with stacked spectra for a single isotope

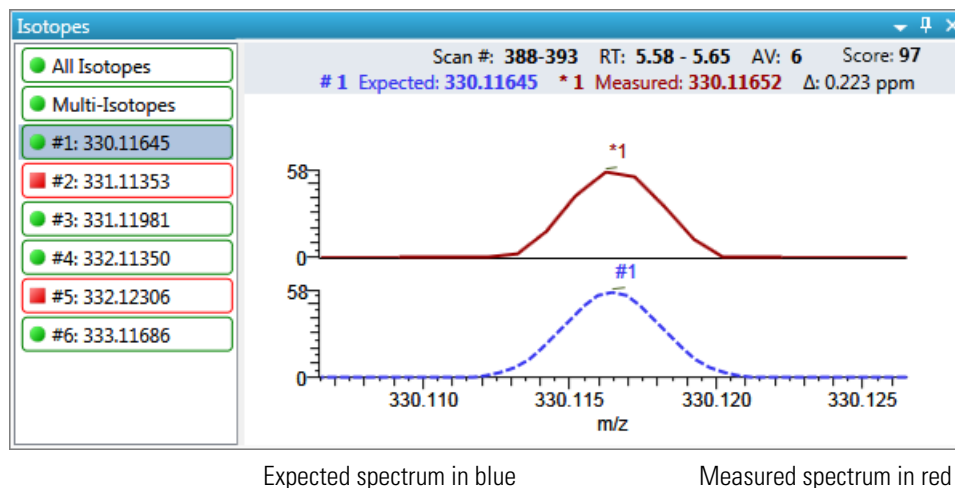
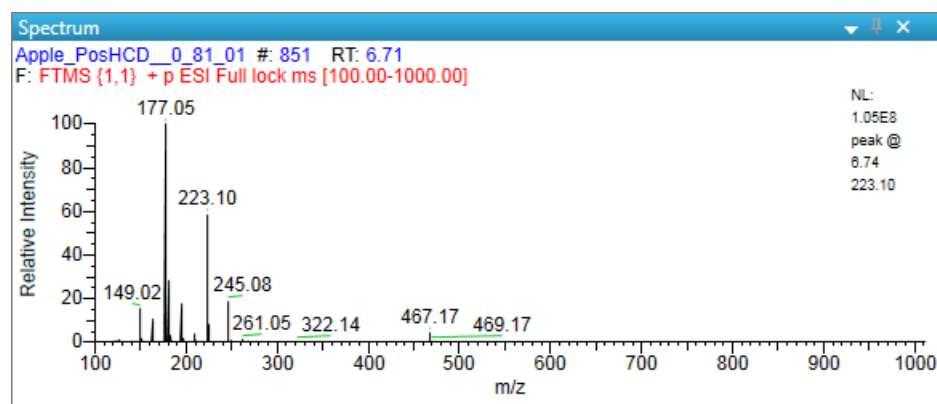


Table 76. Isotopes pane shortcut menu commands

Command	Description
Reset Scaling	Resets the original scaling after a zoom operation.
Copy to Clipboard	Copies the graphic display to the Clipboard.
Display Overlay Spectra Display Stack Spectra	Overlays the two spectrum displays, or stacks the simulated spectrum and the peak apex spectrum.
Show/Hide Noise Label	Adds a noise label to each peak. Expected isotope peaks (displayed in blue) do not display a noise label.
Show/Hide Resolution Label	Adds a resolution label to each peak. Expected isotope peaks (displayed in blue) do not display a resolution label.

Spectrum Pane

When data is available, the plot width is the full mass range in the raw data file. Otherwise, the application scales the width to the scan range.

Figure 150. Spectrum pane**Table 77.** Spectrum pane shortcut menu commands

Command	Description
Reset Scaling	Resets the original scaling after a zoom operation.
Copy to Clipboard	Copies the graphic display to the Clipboard.

Working in the Report View for Unknown Screening Batches

The Report View displays example reports for the current batch. You must have an open batch to use the features in the Report View.

Follow these procedures:

- [To open the Report View](#)
- [To preview a report](#)
- [To generate a report as a PDF, an Excel, or a CSV file](#)
- [To print a report](#)
- [To display a generated report](#)
- [To edit a report template](#)
- [To create a new report template](#)

❖ To open the Report View

Click **Report View** in the navigation pane.



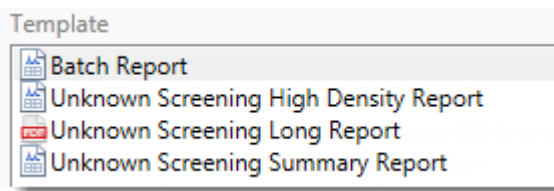
The application opens the [Report View](#).

❖ To preview a report

1. In the Template pane, select a report template.

The template list shows all the report templates that you configured in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.

Figure 151. Example template list for unknown screening batches




2. Click **Preview**, .

The application opens the Report Designer, showing the report information for the current batch in the selected report template format.

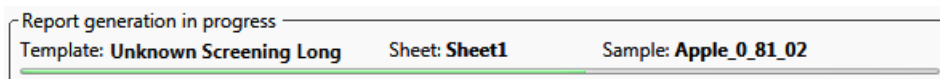
For details about using the Report Designer, see [Chapter 7, “Using the Report Designer.”](#)

❖ **To generate a report as a PDF, an Excel, or a CSV file**

1. In the Template pane, select a report template.
2. Select the check box for each of the file types that you want to create: **PDF**, **Excel**, or **CSV**.
3. Click **Generate**,  **Generate**.


The application does the following:

- Displays a green progress bar as it generates the reports.



- Creates a report for the current batch as a PDF, an Excel, or a CSV file, using the selected report template format.
- Adds information about the generated report to the Generated Reports pane.
For details about the Generated Reports pane, see [Report View](#).
- Saves the report files to the ...\\TraceFinderData\\4.0\\Projects*batch*\\ReportOutput folder.

❖ **To print a report**

1. In the Template pane, select a report template.
2. Select the check box for the **Print** file format.
3. Click **Generate**,  **Generate**.

The application does the following:

- Creates a report for the current batch using the selected report template format.
- Prints the report to your default printer.
- Adds information about the generated report to the Generated Reports pane.
For details about the Generated Reports pane, see [Report View](#).
- Saves the report files to the ...\\TraceFinderData\\4.0\\Projects*batch*\\ReportOutput folder.

❖ **To display a generated report**

In the Generated Reports pane, click **View** for the report that you want to see.

Figure 152. Generated Reports pane showing a PDF report

Generated Reports				
Template	Rule	Sample	Output	Generated Report File
Batch Report	Batch		pdf	View... Batch Report_20140113092939.pdf (3 pages)

Opens the generated file.

The application opens the output file.

❖ To edit a report template

Note You cannot edit report templates that were provided with the TraceFinder application; however, you can open a TraceFinder template, make changes, and save it to a new template name.

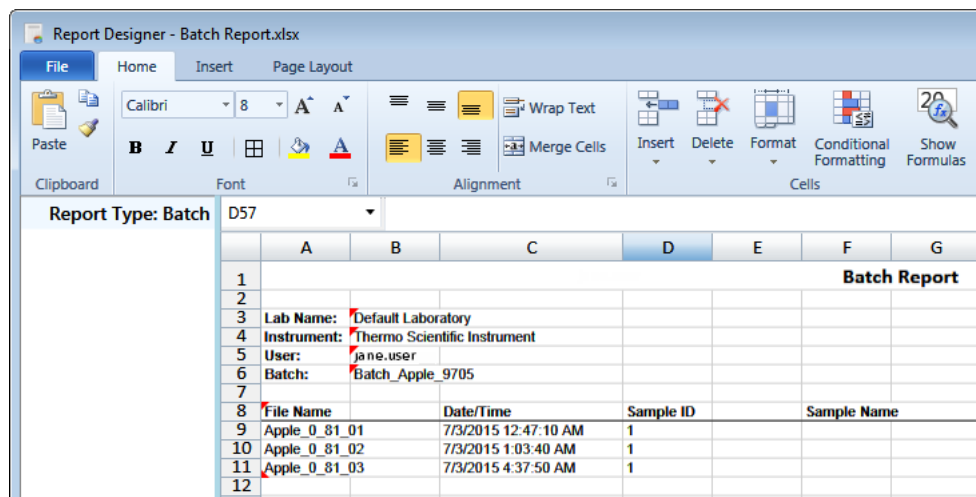
1. In the Template pane, select a report template.

2. Click **Open**, .

The application opens the Report Designer showing the template in an Excel spreadsheet. See [Report Designer showing the template for the selected report](#).

Note When user security is activated, you must have Template Editing permission to edit report templates created by your laboratory. If the Open button is not active, user security is activated and you do not have Template Editing permission.

Figure 153. Report Designer showing the template for the selected report



3. Use the features in the Report Designer to edit the template.

See [Chapter 7, “Using the Report Designer.”](#)

4. When you finish your changes, choose **File > Save** from the Report Designer menu bar.

❖ To create a new report template

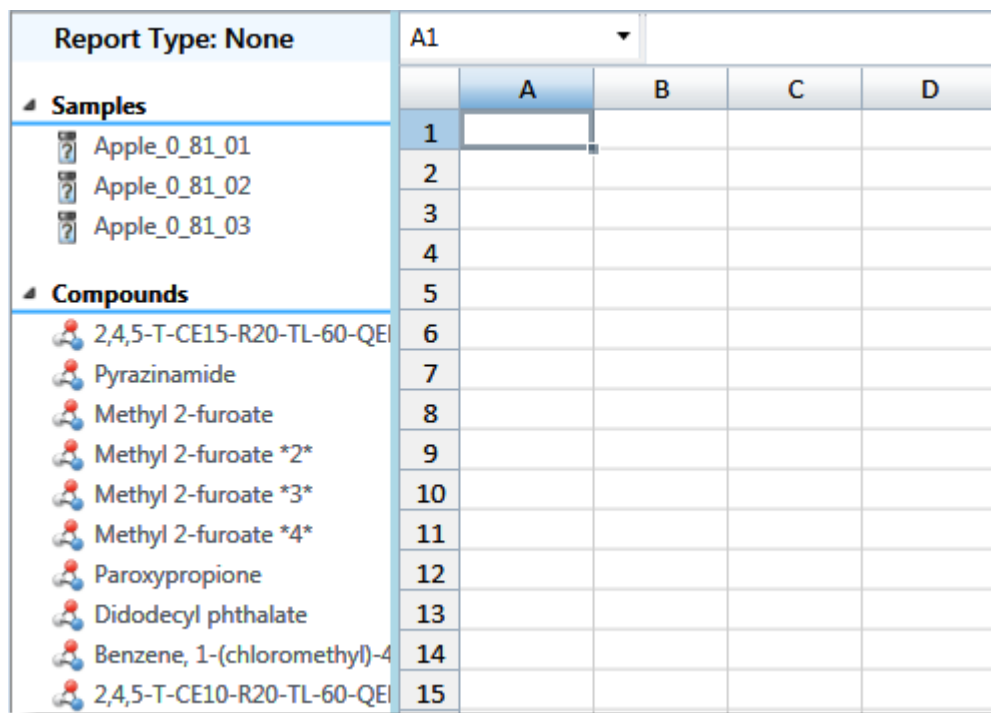
1. Click **New**, .

The application opens the Report Designer showing an empty template in an Excel spreadsheet. See [Report Designer showing a new, empty template](#).

The Report Type is None.

In the left pane, the spreadsheet lists all samples in the current batch and all compounds in the method used for the batch.

Figure 154. Report Designer showing a new, empty template



Report Type: None		A1			
		A	B	C	D
1					
2					
3					
4					
5					
6					
7					
8					
9					
10					
11					
12					
13					
14					
15					

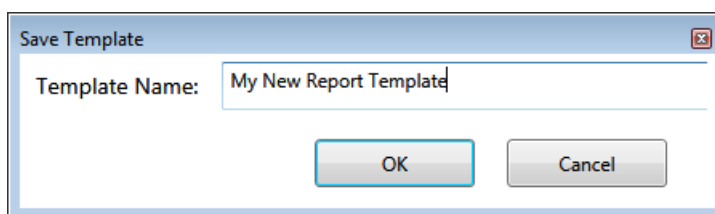
2. Use the features in the Report Designer to create the report template.

See [Chapter 7, “Using the Report Designer.”](#)

3. When you finish your changes, choose **File > Save** from the Report Designer menu bar.

The Save Template dialog box opens.

Figure 155. Save Template dialog box



Save Template

Template Name: My New Report Template

OK Cancel

4. Type a name for the new report template and click **OK**.

Report View

Use the features in the Report View to display example reports for the current batch.

Figure 156. Report View

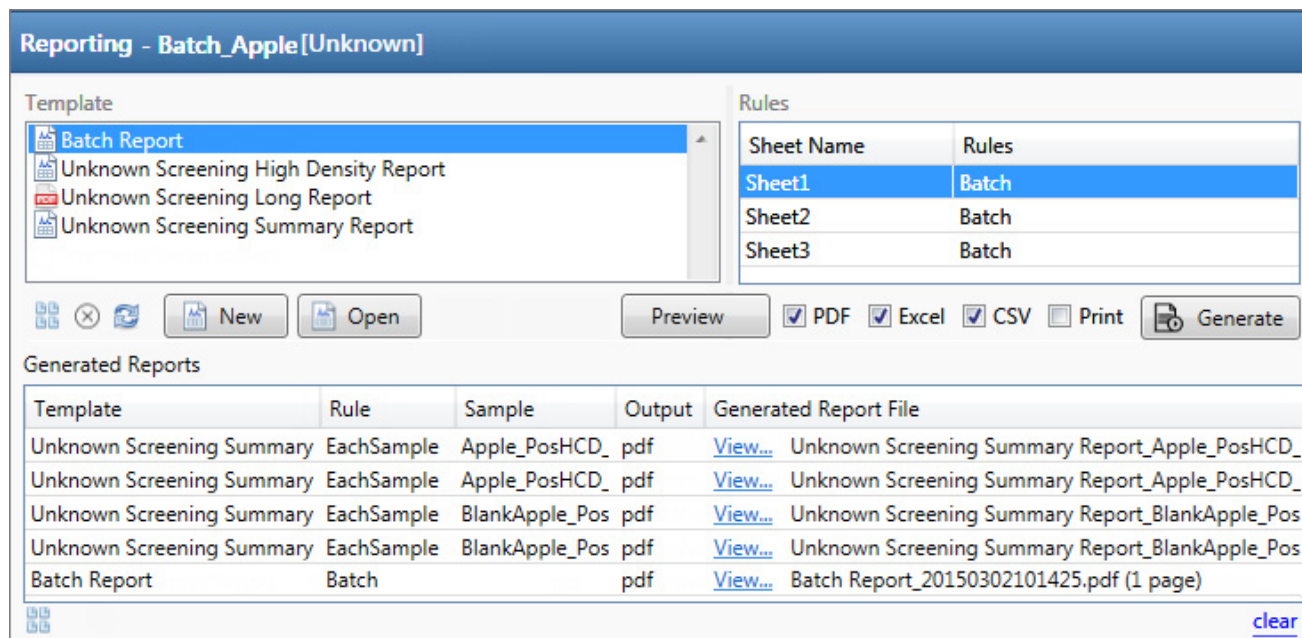





Table 78. Report View parameters (Sheet 1 of 2)

Parameter	Description
Template	
	Displays all available unknown screening report templates.
Rules	
Sheet Name	Specifies each sheet in the report.
Rules	Specifies the type of data used in each sheet in the selected report. <ul style="list-style-type: none"> • Batch • EachSample • SampleType: <i>SampleType</i> • CompoundType: <i>CompoundType</i> • SampleCustomFormula:
Buttons	
View Report Templates	Displays the C:\TraceFinderData\4.0\Templates\ReportTemplates folder that contains all report templates.
Open	Opens the selected report template in the Report Designer.
New	Opens a blank report template in the Report Designer.

Table 78. Report View parameters (Sheet 2 of 2)

Parameter	Description
 Preview	Opens the Report Designer showing the report information for the current batch in the selected report template format.
PDF	Writes the generated report to a PDF file in the ...\\TraceFinderData\\4.0\\Projects\\ <i>batch</i> \\ReportOutput folder.
Excel	Writes the generated report to a PDF file in the ...\\TraceFinderData\\4.0\\Projects\\ <i>batch</i> \\ReportOutput folder.
CSV	Saves the generated report as a PDF file in the ...\\TraceFinderData\\4.0\\Projects\\ <i>batch</i> \\ReportOutput folder. When the report contains multiple sheets, the application writes each sheet as a separate CSV file.
Print	Prints the generated report to your default printer.
 Generate	Generates the selected type of reports for the current batch using the selected report template.
Generated Reports	
Template	Report template used for the report. See Example template list for unknown screening batches .
Rule	Type of data used in each sheet of the report. See Rules .
Sample	For sample-level reports, the name of each sample in the report.
Output	Type of output specified for the report: PDF, Excel, CSV, or Print.
Generated Report File	Lists the output file name for each report in the ...\\TraceFinderData\\4.0\\Projects folder.
View	Displays the generated output file.
 View Generated Reports	Displays the C:\\TraceFinderData\\4.0\\Projects folder that contains all report outputs.
Clear	Removes all reports from the Generated Reports display. This does not delete the reports from the C:\\TraceFinderData\\4.0\\Projects folder.

Working in the Local Method View for Unknown Screening Batches

A local method is a copy of a master method associated with a batch. You can edit only the local copy of the method, or you can edit the master method and overwrite the local copy with the edited master method.

In the Local Method view, you can edit the local method parameters. A local method is a copy of a master method associated with a batch. Local methods are named *Batch_MasterMethod*.

❖ To open the Local Method View

1. Click **Analysis** in the navigation pane.
2. Click **Local Method**.



The [Local Method View](#) for the currently selected batch opens.

You can edit many of the method parameters in a local method. Editing the local method does not affect parameters in the master method.

For detailed descriptions of method parameters, refer to Chapter 6, “Using the Method Development Mode for Unknown Screening Methods,” in the *TraceFinder Lab Director User Guide*.

3. Enter any local changes to the method.
4. When you have finished editing the local method, choose **File > Save**.
5. To process the batch or create new reports with the edited local method, return to the Batch View and submit the batch.

❖ To overwrite the local method with the master method in the Batch View

In the Batch View, click **Update**.



The application overwrites the local method with the master method of the same name. You can use this feature to overwrite an edited local method with the original master method or to overwrite the local method with an updated master method.

6 Using the Analysis Mode for Unknown Screening Batches

Working in the Local Method View for Unknown Screening Batches

Figure 157. Local Method View

The screenshot displays the 'Local Method View' for 'Apple_batch1'. The interface is divided into a left sidebar and a main content area.

Left Sidebar:

- Analysis** (dropdown menu)
- Batch View** (button)
- Data Review** (button)
- Report View** (button)
- Local Method** (dropdown menu, currently expanded)
 - Acquisition** (button)
 - Unknown Screening** (button)
 - Processing** (button)
 - Peak Detection Settings** (button)
 - Reports** (button)

Main Content Area:

Local Method View - Apple_batch1

Master method: [Method Unknown Screening](#)

Lab Name:

Assay type:

Injection Volume:

Mass Precision:

Instrument Method:

Notes:

Using the Report Designer

Use the features in the Report Designer to create or edit report templates.

Contents

- [Editing a Template](#)
- [Toolbar Reference](#)
- [Plot Display Properties](#)
- [Quick Tips](#)

You can create new report templates or edit your laboratory's report templates. You cannot edit report templates that were provided with the TraceFinder application; however, you can open a TraceFinder template, make changes, and save it to a new template name.

When user security is activated, you must have Template Editing permission to edit report templates created by your laboratory. To use the administration tools to set Reports permissions, refer to the instructions in the *TraceFinder Administrator Console User Guide*.

With the Report Designer, you can do any of the following to create or edit a report template:

- Make changes to the template as you would in an Excel application.
- Add or remove graphics.
- Add repeated areas of many types.
- Format fonts and colors.
- Add or edit formulas.
- Edit template text and headers.

In addition to the procedures in this chapter, you can view video instructions at <http://mytracefinder.com/plugins/reporting/>.

Note To edit or create a template, you must have an open batch.

Editing a Template

Use the Report Designer to add or remove graphics, add tables, add repeated areas, and format headings, data, or cells.

Follow these procedures:

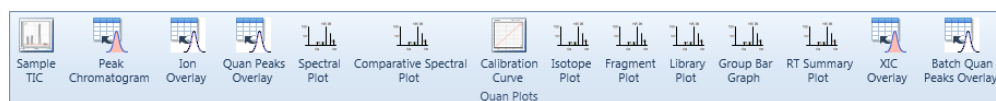
- [To insert graphics into a report template](#)
- [To resize an inserted graphic](#)
- [To move an inserted graphic](#)
- [To insert a table](#)
- [To remove an inserted table](#)
- [To make changes to a table](#)
- [To format header text in the spreadsheet](#)
- [To format data in the report grid](#)
- [To format cells in the report grid](#)
- [To add repeated areas](#)
- [To insert a new sheet in a report](#)

❖ To insert graphics into a report template

1. Click the **Insert** tab.

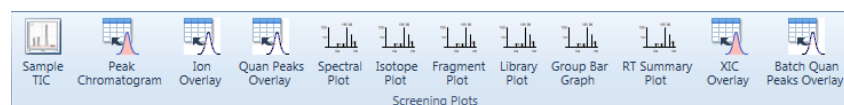
The Insert page displays the available objects that you can insert in a quantitation, a target screening, or an unknown screening report template.

Figure 158. Quan Plots features for quantitation batch report templates



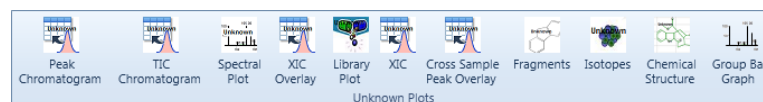
For a description of each graphic element, see [Insert Toolbar – General Features](#).

Figure 159. Screening Plots features for target screening batch report templates



For a description of each graphic element, see [Insert Toolbar – Target Screening Reports](#).

Figure 160. Unknown Plots features for unknown screening batch report templates



For a description of each graphic element, see [Insert Toolbar – Unknown Screening Reports](#).

Note When the method for a quantitation or target screening batch includes unknown screening, the Insert features for the report templates include Unknown Plots.

2. In the template grid, drag to select cells where you want to place the item.

	A	B	C	D
1				
2				
3				
4				
5				

Note You can also select a single row, and the application inserts the upper left corner of the graphic at the first cell, using a minimum default size.

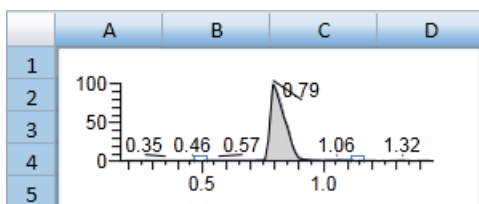
3. Select a sample from the list of samples in the left pane.

The application uses graphical data from the selected sample when you insert a graphical plot.

4. Click a graphical plot in the toolbar to insert into the selected cells.

The application inserts the selected plot in the template grid.

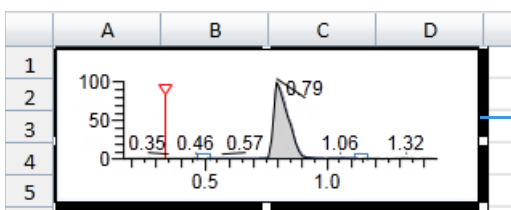
Figure 161. Inserted peak chromatogram



❖ To resize an inserted graphic

1. Right-click the graphic to make it editable.

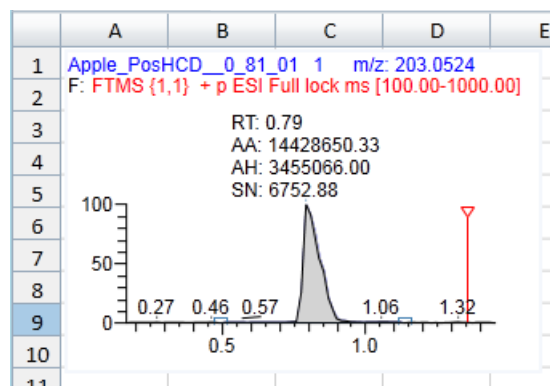
Figure 162. Editable graphic



The border indicates that the graphic is editable.

- Grab a corner of the graphic and stretch it to the new size.

Figure 163. Resized graphic

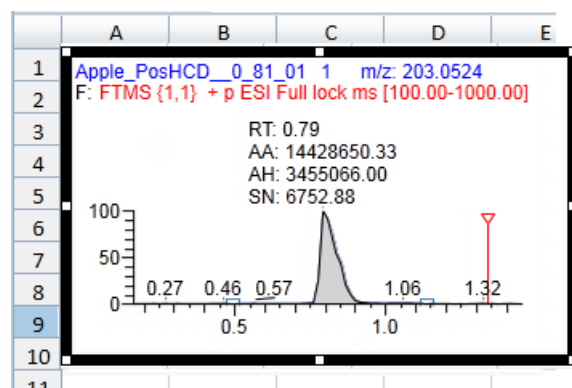


Note In this example, the original graphic was too small to display all the chromatogram data. Stretching the graphic reveals the sample name, mass, filter, and labeling.

❖ To move an inserted graphic

- Right-click the graphic to make it editable.

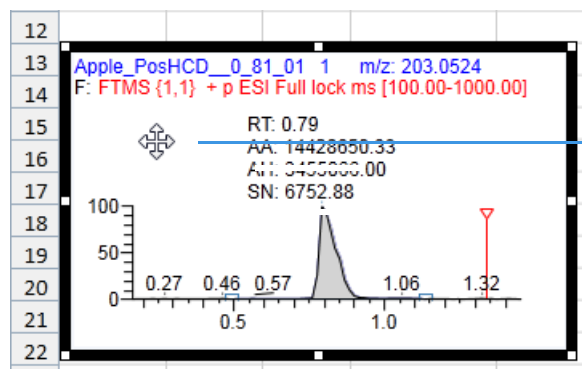
Figure 164. Editable graphic



The border indicates that the graphic is editable.

- Click the graphic (notice the cursor change), and drag it to a new location.

Figure 165. Moved graphic



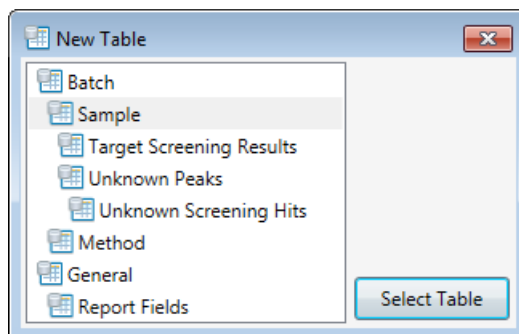
Cursor for moving a graphic

❖ To insert a table

1. Select a row in the report grid where you want to insert the table.
2. Click the **Insert** tab, and then click the **Table** icon,



The New Table dialog box opens.



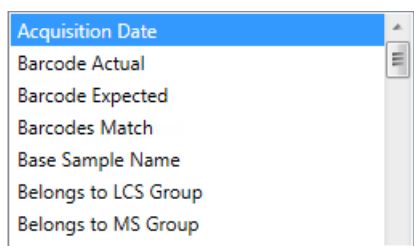
3. Select the type of table that you want to insert and click **Select Table**.

The [Edit Table Dialog Box](#) opens.

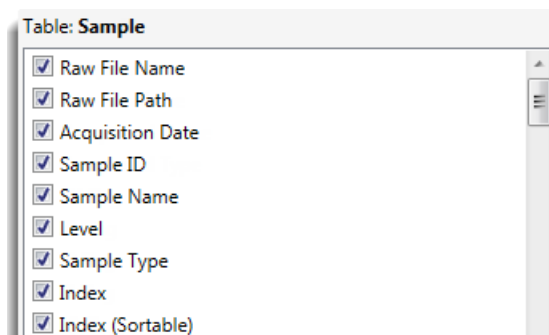
The dropdown list in the upper left corner displays the data type for the table you selected.



The left pane displays an alphabetical list of all available fields for the selected data type that you can include in the table. To display fields for a different data type, see [step 4](#).

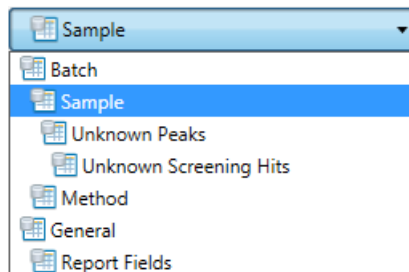


The right pane lists the fields that are included in the new table.




You can select as many additional fields as you want. The table you create displays each selected field as a table column. By default, all available fields for the original data type are included in the table. The first field in the list is the leftmost column in the table. The last field in the list is the rightmost column in the table. To change the order of the columns, see [step 6](#).



4. To add fields to the table from a different data type, do the following:
 - a. Click the dropdown list and select a different data type.



The left pane displays an alphabetical list of all available fields for the new data type.

- b. Select a field in the left pane and then click .

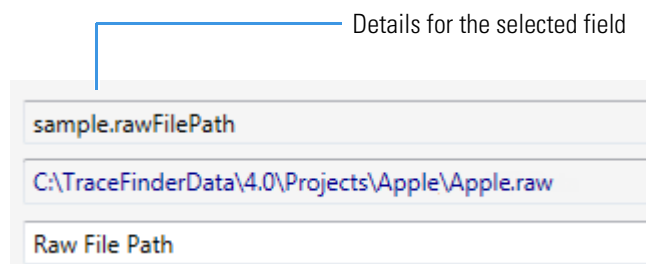
The application moves the selected field to the rightmost pane to indicate that it is included in the inserted table.

5. To select a field that you do not want in the table, in the rightmost pane, clear the check box next to the field name.
6. To move a column up or down in the table, select the field name in the rightmost pane and click the **Up** arrow, , or **Down** arrow, .
7. To edit the header for a table column, select the field in the list on the right, and then type new header text in the Header box in the Field Details area.

The default header name is the same as the field name.

8. To find the formula for a field, select the field.

The application displays the formula in the Formula box.



This video demonstrates how to use the Report Designer to create your own formula:
[Working with Formulas](#).

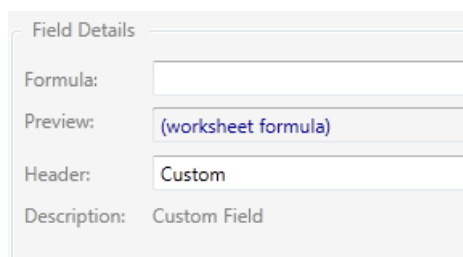
9. To create a custom column, do the following:

- a. Click **Custom**, .

The application adds a Custom field to the fields list.

- b. Select the Custom field in the list.

The application display the details for the new, custom column.



The 'Field Details' dialog box contains the following fields:

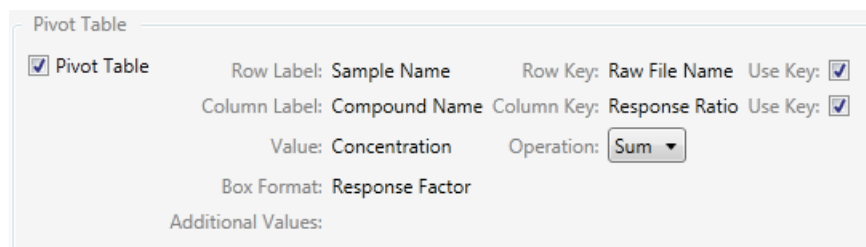
- Formula:** (empty text box)
- Preview:** (worksheet formula)
- Header:** Custom
- Description:** Custom Field

- c. Enter values for the Formula and Header.

10. To create a pivot table, do the following:

- a. Select the **Pivot Table** check box.

The Pivot Table area expands to display the pivot table options.



The 'Pivot Table' dialog box contains the following options:

- ☒ **Pivot Table**
- Row Label: Sample Name
- Row Key: Raw File Name Use Key: ☒
- Column Label: Compound Name
- Column Key: Response Ratio Use Key: ☒
- Value: Concentration
- Operation: Sum
- Box Format: Response Factor
- Additional Values:

- b. Select the **Use Key** check box for the row label.
- c. Select the **Use Key** check box for the column label.
- d. Select the operation that you want to use to calculate the aggregate value: **Sum**, **Count**, **Average**, **First**, or **Last**.

11. When you have made all your table selections, click **OK**.

The application inserts the table into the report template.

❖ To remove an inserted table

1. Select all rows and cells in the table.
2. Right-click and choose **Delete – Delete Cells and Shift Cells Left**.

The application removes the entire table from the report template.

❖ **To make changes to a table**

1. In the table that you want to edit, right-click and choose **Data Table**.

The **Edit Table Dialog Box** opens, displaying all of the available fields for the selected template. The Edit Table dialog box is virtually identical to the dialog box that you used to create a table.

2. Edit the fields for the table and click **OK**.

❖ **To format header text in the spreadsheet**

1. In the Report Designer window, select the header in the spreadsheet table.

To change all headers, select the entire header row.

To change a single header, select the specific header cell.

Tip You can also use the SHIFT key to select sequential cells or the CTRL key to select nonsequential cells anywhere in the grid.

2. Click the **Home** tab.
3. Use the Font icons to change the font format.



❖ **To format data in the report grid**

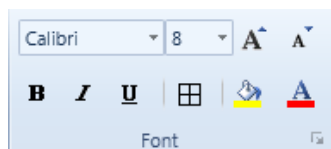
1. Select the data that you want to format.

To format all rows in a column, select the header row.

To format a single cell of data, select only that cell.

Tip You can also use the SHIFT key to select sequential cells or the CTRL key to select nonsequential cells anywhere in the grid.

2. Click the **Home** tab and use the toolbar icons to edit the font or cells as appropriate:



- Change the font or font size.
- Make the selected text bold.
- Make the selected text italics.

- Underline the selected text.
- Apply borders to the currently selected cells.
- Increase or decrease font size.
- Apply color to the background for selected cells.
- Change the font color.

❖ **To format cells in the report grid**

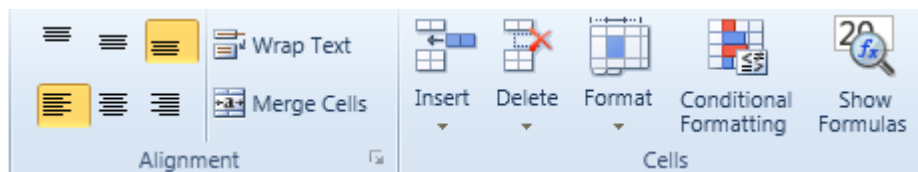
1. Select the cells that you want to modify:

To change all cells in a row, select the entire row.

To change a single cell, select only that cell.

Tip You can also use the SHIFT key to select sequential cells or the CTRL key to select nonsequential cells.

2. Click the **Home** tab and use the Alignment or Cells toolbar icons to edit the cells.



3. To align the cell text to the top, center, or bottom of the cell, click one of these icons:



4. To align the cell text to the left, center, or right of the cell, click one of these icons:



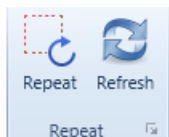
5. To make all contents visible within a cell, click **Wrap Text**.
6. To join selected cells into one cell, click **Merge Cells**.
7. To insert cells, rows, or columns into the template, click **Insert**.
8. To delete rows or columns from the template, click **Delete**.
9. To change the row height or column width, organize sheets, or protect or hide cells, click **Format**.
10. To highlight or emphasize useful cells based on specific criteria, click **Conditional Formatting** to use data bars, color scales, or icon sets.
11. To show formulas for selected cells instead of the resulting value, click **Show Formulas**.

❖ To add repeated areas

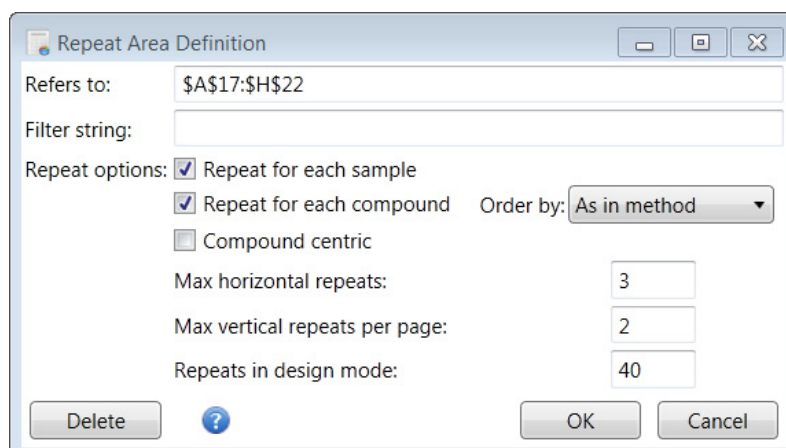
1. To define a repeat area, select the cells that you want to repeat.

You can repeat only one area per worksheet, but you can repeat it many times. You cannot insert data tables beneath a repeated area. See [Example of repeated areas](#).

2. Click the **Insert** tab and then click **Repeat**.



The Repeat Area Definition dialog box opens.

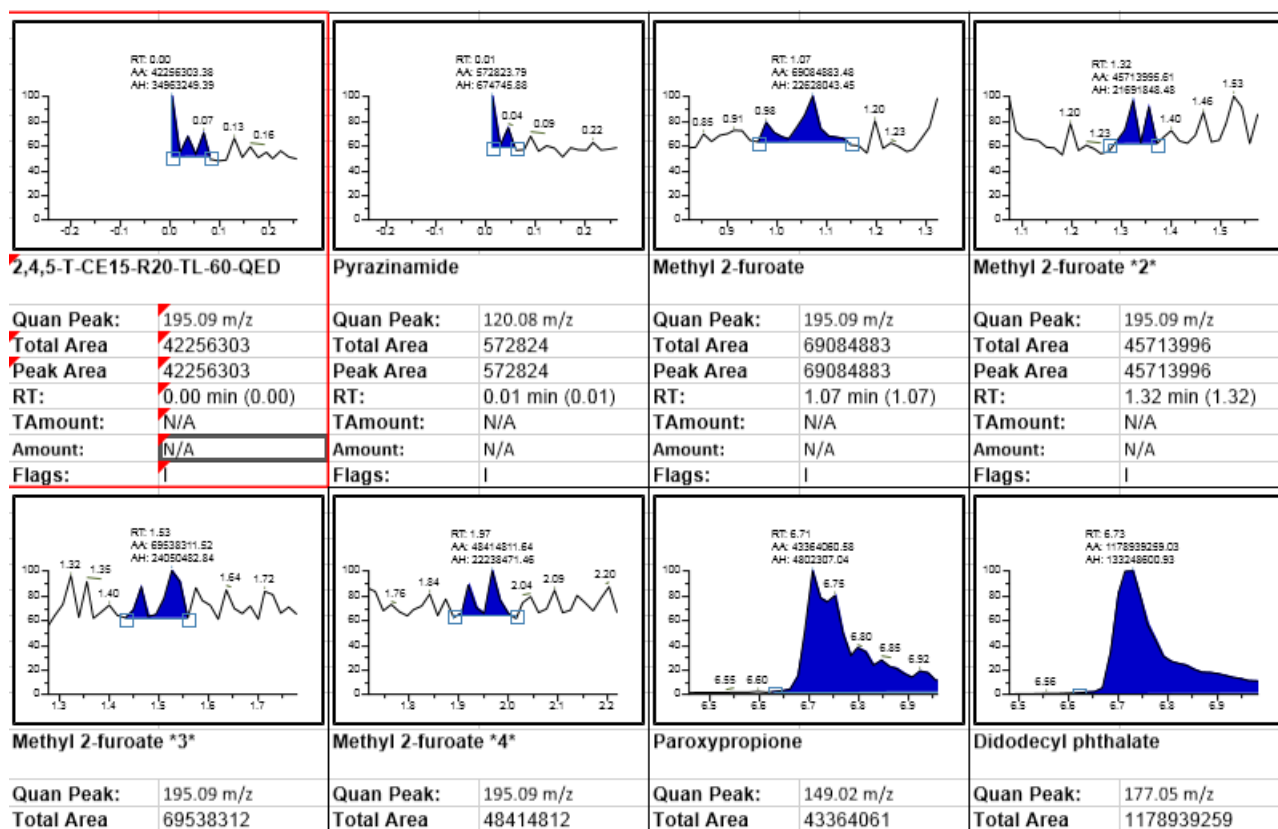


3. To filter the repeat area according to the filter criteria, type the string syntax in the Filter String box.
4. Select to repeat for each sample, for each compound, or for both.
5. Define whether the repeat area repeats first for samples or first for compounds:
 - To have the repeat area repeat first for samples and then for compounds, clear the **Compound Centric** check box.
 - To have the repeat area repeat first for compounds and then for samples, select the **Compound Centric** check box.
6. To define the number of times to repeat horizontally before wrapping to the next row, type a number in the Max Horizontal Repeats box.
7. To define the number of times to repeat vertically before inserting an auto-page break, type a number in the Max Vertical Repeats Per Page box.
8. To define the maximum suggested number of repeats, type a number in the Repeats in Design Mode box.

This maximum number of repeats is intended to limit the size of the report while you are editing the report in the Report Designer. When generating an actual report, the application does not enforce this limit.

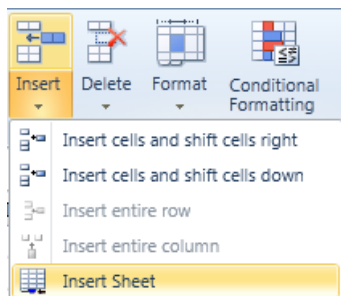
In this example, the repeated area (outlined in red) displays the peak chromatogram, target peak, total area, peak area, retention time, theoretical amount, actual amount, and reported flags, if any, for each compound.

Figure 166. Example of repeated areas



❖ To insert a new sheet in a report

1. Click the **Home** tab.
2. In the Cells area, choose **Insert Sheet** from the Insert menu.



The application adds a new sheet to the report.

Edit Table Dialog Box

Use the options in the Edit Table dialog box to edit the fields in a table.


Figure 167. Edit Table dialog box

The screenshot shows the 'Edit Table' dialog box. At the top, there's a 'Sample' dropdown menu. Below it is a list of fields: Acquisition Date, Base Sample Name, Channel, Comment, Index, Index (Sortable), Injection Volume, Instrument Method, Instrument Method Acquired, Raw File Name (highlighted), and Raw File Path. To the right of this list are navigation buttons: '>', '>>', '<<', '<', '<<<', '>>>', and 'Custom'. Further right is a 'Table: Sample' list with checkboxes for: Raw File Name, Raw File Path, Acquisition Date, Sample ID, Sample Name, Index, Index (Sortable), Vial Position, Injection Volume, Comment, State, Channel, and Base Sample Name. Below these is the 'Field Details' section with 'Formula:' (sample.rawFilename), 'Preview:' (Blank_PosHCD_02), and 'Description:' (The user-defined name of the rawfile.). To the right of this is a 'With Extension' checkbox. Below that is the 'Table Filter' section with 'Table Filter:' (result.compoundType <> "Internal Standard"), 'Preview:', 'Sort By: Sample ID' (with a sort icon), 'Then By:' (with a sort icon), and a 'Repeat Header Per Group' checkbox. At the bottom is the 'Pivot Table' section with a 'Pivot Table' checkbox, 'Row Label: Sample Name', 'Row Key: Raw File Name' (with 'Use Key: [checked]'), 'Column Label: Compound Name', 'Column Key: Response Ratio' (with 'Use Key: [checked]'), 'Value: Concentration', 'Operation: Sum' (in a dropdown), 'Box Format: Response Factor', and 'Additional Values:'. An 'OK' button is at the bottom right.

Table 79. Edit Table dialog box parameters (Sheet 1 of 2)

Parameter	Description
Dropdown List	Specifies the type of table you are editing.
Field List	Lists all available table fields.
Table	Displays the selected table type and the fields included in the table.

Table 79. Edit Table dialog box parameters (Sheet 2 of 2)

Parameter	Description
Field Details	
Formula	Specifies the functions and arguments used to define the field.
Preview	Specifies the name of the raw data file used for the report.
Header	Specifies the header name for the column in the spreadsheet.
Description	Specifies the type of field.
Table Filter	
Table Filter	Specifies the filter used for the sample compounds.
Preview	—
Sort By	Sorts the data by ascending or descending values.
Pivot Table	
Pivot Table	Enables the Pivot Table features.
Row Label	—
Row Key	—
Column Label	—
Column Key	—
Value	—
Operation	Specifies the operation used to calculate the aggregate value. Valid values: Sum, Count, Average, First, Last
Box Format	—
Additional Values	—
 Delete	Deletes the selected table from the template.

Toolbar Reference

The Report Designer includes the following toolbars.

- [Home Toolbar](#)
- [Insert Toolbar – General Features](#)
- [Insert Toolbar – Quantitation Reports](#)
- [Insert Toolbar – Target Screening Reports](#)
- [Insert Toolbar – Unknown Screening Reports](#)
- [Page Layout Toolbar](#)

Home Toolbar

Use the options in the Home toolbar to modify fonts, align cell data, and format cells.

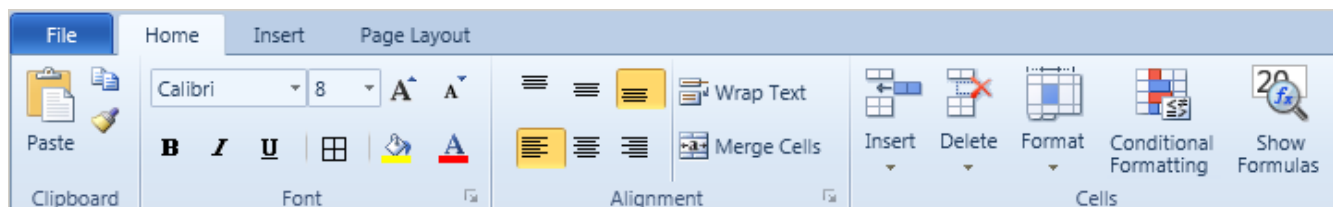


Table 80. Home toolbar options (Sheet 1 of 2)

Option	Description
Clipboard	
Paste	Paste the contents of the Clipboard. You can also paste only formatting or only a formula.
Copy	Copy text or graphics to use in another place.
Font	
Font and size lists	Use the dropdown lists to change the font or font size.
Bold	Make the selected text bold.
Italic	Make the selected text italics.
Underline	Underline the selected text.
Border	Apply borders to the currently selected cells.
Font size	Increase or decrease the font size.
Fill color	Apply color to the background for selected cells.
Font color	Change the font color.
Alignment	
	Align the cell text to the top, center, or bottom of the cell.
	Align the cell text to the left, center, or right of the cell.
Wrap text	Make all contents visible within a cell.

Table 80. Home toolbar options (Sheet 2 of 2)

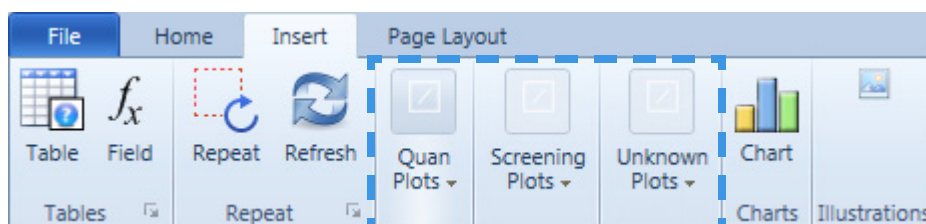
Option	Description
Merge cells	Join selected cells into one cell.
Cells	
Insert	Insert cells, rows, columns, or sheets into a template.
Delete	Delete rows or columns from a template.
Format	Change the row height or column width, organize sheets, or protect or hide cells.
Conditional Formatting	Based on specific criteria, highlight or emphasize useful cells using data bars, color scales, and icon sets.
Show Formulas	Show formulas for selected cells instead of the resulting value.

Insert Toolbar – General Features

Use the features on the Insert page to create and edit tables, formulas, graphic plots, charts, and illustrations.

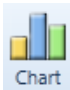
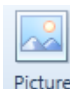
Graphic plots can represent data from quantitative, target screening, or unknown screening data. The Insert page displays the available data plots for the type of batch: quantitation, quantitation plus unknown screening, target screening, target screening plus unknown screening, or unknown screening only.

The following graphic outlines all three plot types for illustrative purposes only.

**Table 81.** Insert toolbar options (Sheet 1 of 2)

Option	Description
Tables	
Table	Create a table to manage and analyze data.
Field	Edit the formula for a field by choosing functions and editing arguments.
Repeat	
Repeat	Repeat text, cell, or graphic elements.
Refresh	Update the view to reflect recent changes.
Quan Plots	Can be stand-alone or used with Unknown plots. For descriptions of quantitation plots, see Insert Toolbar – Quantitation Reports .
Screening Plots	Can be stand-alone or used with Unknown plots. For descriptions of target screening plots, see Insert Toolbar – Target Screening Reports .

Table 81. Insert toolbar options (Sheet 2 of 2)

Option	Description
Unknown Plots	Can be stand-alone or used with either Quan plots or Screening plots. For descriptions of unknown screening plots, see Insert Toolbar – Unknown Screening Reports .
Charts	
	Adds a custom chart. To define the chart, double-click the inserted blank chart.
Illustrations	
	Insert a picture from a file.

Insert Toolbar – Quantitation Reports

Use the options in the Insert toolbar to add graphics, plots, and other objects to a template or report. You can also set up repeating objects and define functions.

For each inserted plot, you can customize the display properties. See [Quantitation and Target Screening Plots](#).

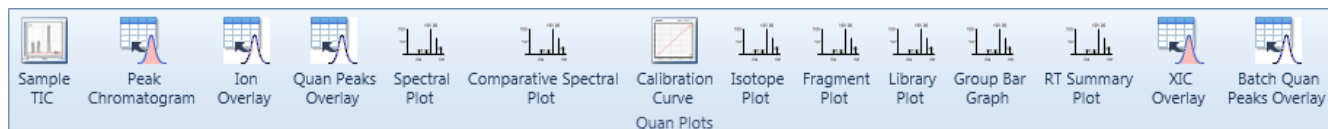


Table 82. Insert toolbar Quantitation options (Sheet 1 of 2)

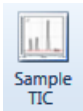
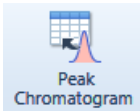
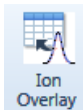

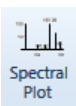

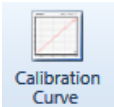





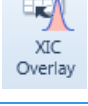
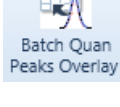
Option	Description
	Adds a sample TIC. For information about sample TIC data in the Data Review results, see Peaks Pane .
	Adds a peak chromatogram. For information about peak chromatogram data in the Data Review results, see Quan Peak .
	Adds an ion overlay. For information about ion overlay data in the Data Review results, see Ion Overlay .
	Adds a quantitation peaks overlay plot. For information about quantitation peaks overlay data in the Data Review results, see Quan Peaks Overlay .

Table 82. Insert toolbar Quantitation options (Sheet 2 of 2)

Option	Description
 <p>Spectral Plot</p>	<p>Adds a spectral plot.</p> <p>For information about spectral data in the Data Review results, see Spectra.</p>
 <p>Comparative Spectral Plot</p>	<p>Adds a comparative spectral plot.</p> <p>For information about comparative spectral data in the Data Review results, see Spectra.</p>
 <p>Calibration Curve</p>	<p>Adds a calibration curve plot.</p> <p>For information about calibration curve data in the Data Review results, see Calibration Curve.</p>
 <p>Isotope Plot</p>	<p>Adds an isotope plot to display the number of isotopes found, the score, a pass/fail flag, and a plot of the isotopes.</p> <p>For information about isotope plot data in the Data Review results, see Isotope.</p>
 <p>Fragment Plot</p>	<p>Adds a fragment plot.</p> <p>For information about fragment plot data in the Data Review results, see Fragments.</p>
 <p>Library Plot</p>	<p>Adds a library plot.</p> <p>For information about library plot data in the Data Review results, see Library Match.</p>
 <p>Group Bar Graph</p>	<p>Adds a group averages plot.</p> <p>For information about group averages data in the Data Review results, see Group Averages.</p>
 <p>RT Summary Plot</p>	<p>Adds an RT summary plot.</p> <p>For information about retention time summary data in the Data Review results, see Retention Time Summary.</p>
 <p>XIC Overlay</p>	<p>Adds an XIC summary plot.</p> <p>For information about XIC summary data in the Data Review results, see XIC Overlay.</p>
 <p>Batch Quan Peaks Overlay</p>	<p>Adds a Batch Quan Peaks Overlay plot that displays all quantitative peaks in the batch.</p>

Insert Toolbar – Target Screening Reports

Use the options in the Insert toolbar to add graphics, plots, and other objects to a template or report. You can also set up repeating objects and define functions.

For each inserted plot, you can customize the display properties. See [Quantitation and Target Screening Plots](#).

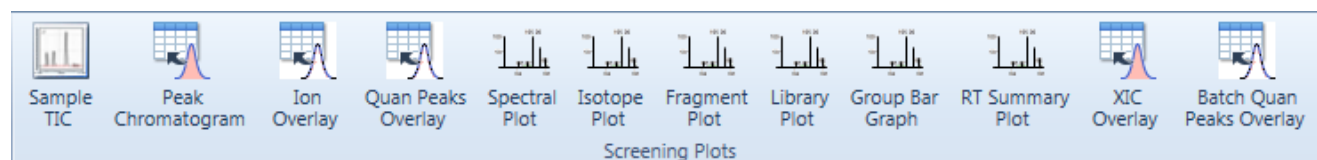


Table 83. Insert toolbar Target Screening options (Sheet 1 of 2)

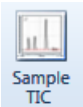
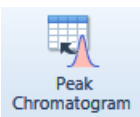
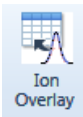
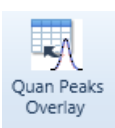

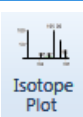
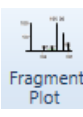



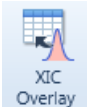
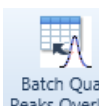
Option	Description
	Adds a sample TIC plot.
	Adds a peak chromatogram plot. For information about peak chromatogram data in the Data Review results, see Chromatogram Pane .
	Adds an ion overlay plot.
	Adds a quantitation peaks overlay plot.
	Adds a spectral plot. For information about spectral plot data in the Data Review results, see Spectrum .
	Adds an isotope plot to display the number of isotopes found, the score, a pass/fail flag, and a plot of the isotopes. For information about isotope plot data in the Data Review results, see Isotopes .
	Adds a fragment plot. For information about fragment plot data in the Data Review results, see Fragments .
	Adds a library plot. For information about library plot data in the Data Review results, see Library .

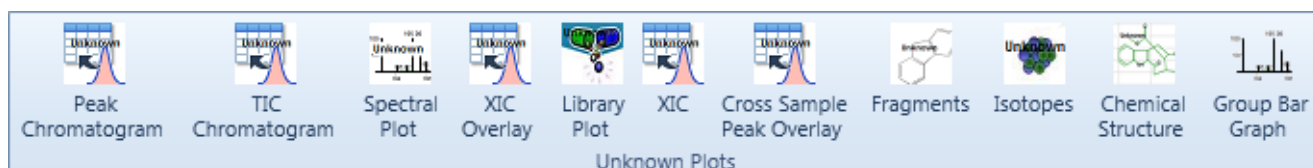
Table 83. Insert toolbar Target Screening options (Sheet 2 of 2)

Option	Description
	Adds a group averages plot. For information about group averages data in the Data Review results, see Groups .
	Adds an RT summary plot.
	Adds an XIC summary plot.
	Adds a Batch Quan Peaks Overlay plot that displays all quantitative peaks in the batch.

Insert Toolbar – Unknown Screening Reports

Use the options in the Insert toolbar to add graphics, plots, and other objects to a template or report. You can also set up repeating objects and define functions.

For each inserted plot, you can customize the display properties. See [Unknown Screening Plots](#).

**Table 84.** Insert toolbar Unknown Screening options (Sheet 1 of 2)

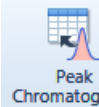
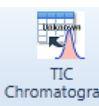

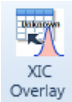
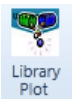
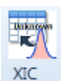
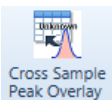
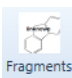

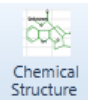

Option	Description
	Adds a peak chromatogram. For information about peak chromatogram data in the Data Review results, see Peak Chromatogram Pane .
	Adds a TIC chromatogram. For information about TIC chromatogram data in the Data Review results, see TIC Chromatogram Pane .
	Adds a spectral plot. For information about spectral plot data in the Data Review results, see Spectrum Pane .

Table 84. Insert toolbar Unknown Screening options (Sheet 2 of 2)

Option	Description
	<p>Adds an XIC overlay plot.</p> <p>For information about XIC overlay data in the Data Review results, see XIC Overlay Pane.</p>
	<p>Adds a library plot.</p> <p>For information about library plot data in the Data Review results, see Library Search Pane.</p>
	<p>Adds an XIC plot.</p> <p>For information about XIC plot data in the Data Review results, see XIC Pane.</p>
	<p>Adds a cross-sample peak overlay plot.</p> <p>For information about cross-sample peak overlay data in the Data Review results, see Cross Sample Peak Overlay Pane.</p>
	<p>Adds a fragments plot.</p> <p>For information about fragments plot data in the Data Review results, see Fragments Pane.</p>
	<p>Adds an isotopes plot.</p> <p>For information about isotopes plot data in the Data Review results, see Isotopes Pane.</p>
	<p>Adds a chemical structure plot.</p> <p>For information about chemical structure plot data in the Data Review results, see Chemical Structure Pane.</p>
	<p>Adds a group averages plot.</p> <p>For information about group averages data in the Data Review results, see Group Averages Pane.</p>

Page Layout Toolbar

Use the options in the Page Layout toolbar to adjust margins, orientation, paper size, and page breaks. To see your changes, click **Print Preview**.

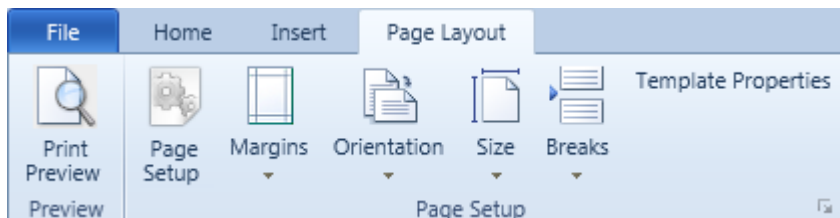
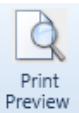
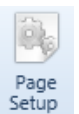

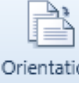

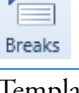


Table 85. Page Layout toolbar options

Option	Description
 Print Preview	View the report before printing it.
 Page Setup	Define page details.
 Margins	Select page margins for the current view or the entire document.
 Orientation	Switch the pages from portrait to landscape view.
 Size	Choose a page size for the current view or the entire document.
 Breaks	Specify where a new page begins in the printed copy.
Template Properties	<p>Open the Template Properties dialog box where you can specify the expected batch type for the report template.</p> <p>The Template Properties dialog box also displays the Used Range so that you can identify when there is empty white space in the report template. Extra white space to the right or left of the report content can cause unwanted page breaks.</p>

Plot Display Properties

For plot types that you insert in the spreadsheet, you can change the display properties.

Note This topic shows examples of all plot types; however, the following plots do not have editable plot properties: Calibration Curve, Group Bar Graph, and RT Summary.

❖ To change the plot display properties

1. Insert a plot in the spreadsheet.
2. Right-click the plot to select it.
3. Right-click the plot again to display the shortcut menu.
4. Choose **Properties**.

The Plot Properties dialog box opens, displaying the editable display properties for the selected plot type.

The following topics show the plots in the report templates and describe their properties.

- [Quantitation and Target Screening Plots](#)
- [Unknown Screening Plots](#)

Quantitation and Target Screening Plots

To customize the display in the quantitation and target screening reports, use the Plot Properties dialog box for each plot.

- [Sample TIC Plot](#)
- [Peak Chromatogram Plot](#)
- [Ion Overlay Plot](#)
- [Quan Peaks Overlay Plot](#)
- [Spectral Plot](#)
- [Comparative Spectral Plot \(quantitation reports only\)](#)
- [Calibration Curve Plot \(quantitation reports only\)](#)
- [Isotope Plot, Fragment Plot, or Library Plot](#)
- [Group Bar Graph Plot](#)
- [RT Summary Plot](#)
- [XIC Overlay Plot](#)
- [Batch Quan Peaks Overlay Plot](#)

Sample TIC Plot

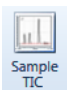
Use the  icon to insert a sample TIC Chromatogram plot.

Figure 168. TIC Chromatogram plot

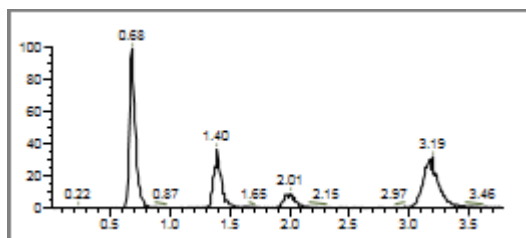


Figure 169. Plot Properties dialog box for a TIC Chromatogram plot

Table 86. Plot Properties parameters for a TIC Chromatogram (Sheet 1 of 2)

Command	Description
TIC Polarity	Specifies the polarity of adducts displayed in the plot. All: Displays positive, negative, and analog adducts. Positive: Displays only positive adducts. Negative: Displays only negative adducts. Analog: Displays only analog adducts.
Y-Axis Type	
Relative	Displays the y-axis scale from 0 through 100.
Absolute	Displays the y-axis scale from 0 to the actual value of the most intense peak in the group.

Table 86. Plot Properties parameters for a TIC Chromatogram (Sheet 2 of 2)

Command	Description
Label in Scientific Notation	Displays the <i>y</i> -axis scale in scientific notation.
Labels	
Show Label for Y-Axis	Displays the relative intensity value on the <i>y</i> axis.
Show Label for X-Axis	Displays the retention time (RT) value on the <i>x</i> axis.
Show Header	Displays the raw data file path at the top of the plot.
Show Grid Lines	Displays the grid lines in the plot background.
Peak Plot Labels	
RT	Displays the retention time (RT) for each peak in the plot.
Show All Apex Times	Displays all surrounding apex times in the plot.
Nominal Level	Displays the nominal level (NL) values for each peak in the plot.

Peak Chromatogram Plot

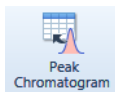
Use the  icon to insert a Peak Chromatogram plot.

Figure 170. Peak Chromatogram plot

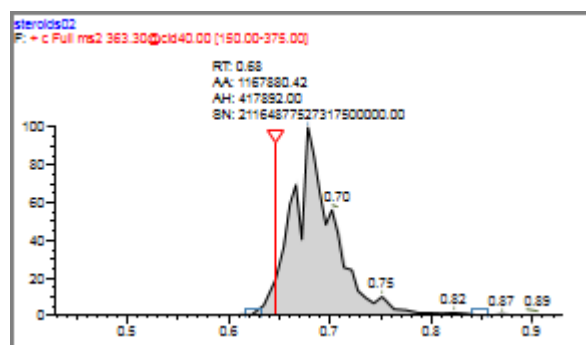
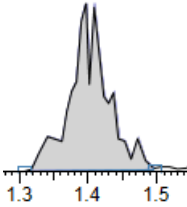
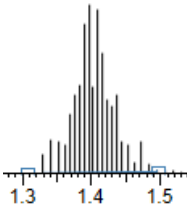


Figure 171. Plot Properties dialog box for Peak Chromatogram plot

Table 87. Plot Properties parameters for a Peak Chromatogram plot (Sheet 1 of 2)

Command	Description
Peak Type	Specifies the type of peak to display in the plot: <ul style="list-style-type: none"> • Quan • Confirming • InternalStandard • InternalStandardConfirming
Integration: Always Show Processed Peak	Specifies that the plot always displays the processed peak.
Quan Peak Index	Valid values: 1 through 9
Time Range	Specifies that the plot display one of the following: <ul style="list-style-type: none"> • View Width: Zooms the display on the peak of interest. • Full Window: Displays the full width of the scan.
Y-Axis Type	
Relative	Displays the y-axis scale from 0 through 100.
Absolute	Displays the y-axis scale from 0 to the actual value of the most intense peak in the group.

Table 87. Plot Properties parameters for a Peak Chromatogram plot (Sheet 2 of 2)

Command	Description
Label in Scientific Notation	Displays the <i>y</i> -axis scale in scientific notation.
Labels	
Show Label for Y-Axis	Displays the relative intensity value on the <i>y</i> axis.
Show Label for X-Axis	Displays the retention time (RT) value on the <i>x</i> axis.
Show Grid Lines	Displays the grid lines in the plot background.
Peak Plot Labels	
RT	Displays the retention time (RT) label in the peak chromatogram plot.
Area	Displays the peak area (AA) label in the peak chromatogram plot.
Height	Displays the peak height (AH) label in the peak chromatogram plot.
Signal to Noise	Displays the signal-to-noise (SN) label in the peak chromatogram plot.
Show All Apex Times	Displays all surrounding apex times in the plot.
Sample Filename	Displays the name of the sample file.
Compound Name	Displays the compound name.
Nominal Level	Displays the nominal level (NL) values for each peak in the plot.
Filter	Displays the filter used for the scan. The plot displays the filter name below the raw data file name.
Plot Style	
Point To Point	Displays chromatograms in the following format: 
Stick	Displays chromatograms in the following format: 

Ion Overlay Plot

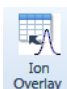
Use the  icon to insert an Ion Overlay plot.

Figure 172. Ion Overlay plot

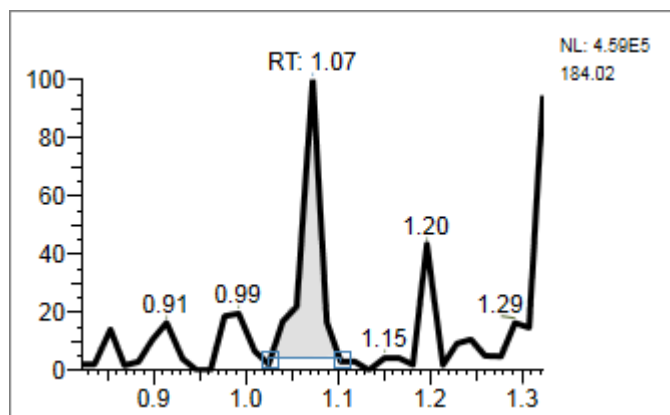


Figure 173. Plot Properties dialog box for Ion Overlay plot

Table 88. Plot Properties parameters for an Ion Overlay plot (Sheet 1 of 2)

Command	Description
Peak Type	Specifies the type of peak to display in the plot: Quan or InternalStandard
Quan Peak Index	Valid values: 1 through 9

Table 88. Plot Properties parameters for an Ion Overlay plot (Sheet 2 of 2)

Command	Description
Time Range	Specifies that the plot display one of the following: <ul style="list-style-type: none"> View Width: Zooms the display on the peak of interest. Full Window: Displays the full width of the scan.
Y-Axis Type	
Relative	Displays the <i>y</i> -axis scale from 0 through 100.
Absolute	Displays the <i>y</i> -axis scale from 0 to the actual value of the most intense peak in the group.
Label in Scientific Notation	Displays the <i>y</i> -axis scale in scientific notation.
Labels	
Show Label for Y-Axis	Displays the relative intensity value on the <i>y</i> axis.
Show Label for X-Axis	Displays the retention time (RT) value on the <i>x</i> axis.
Show Header	Displays the raw data file name and compound name.
Show Grid Lines	Displays the grid lines in the plot background.

Quan Peaks Overlay Plot

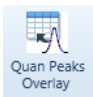
Use the  icon to insert a quantitation peaks overlay plot.

Figure 174. Quan Peaks Overlay plot

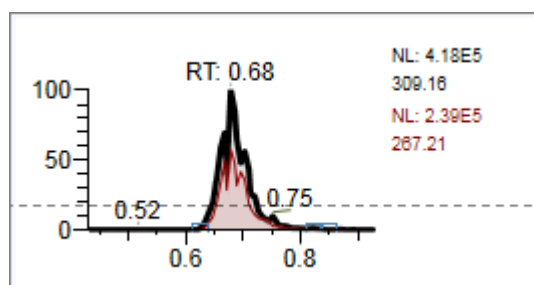
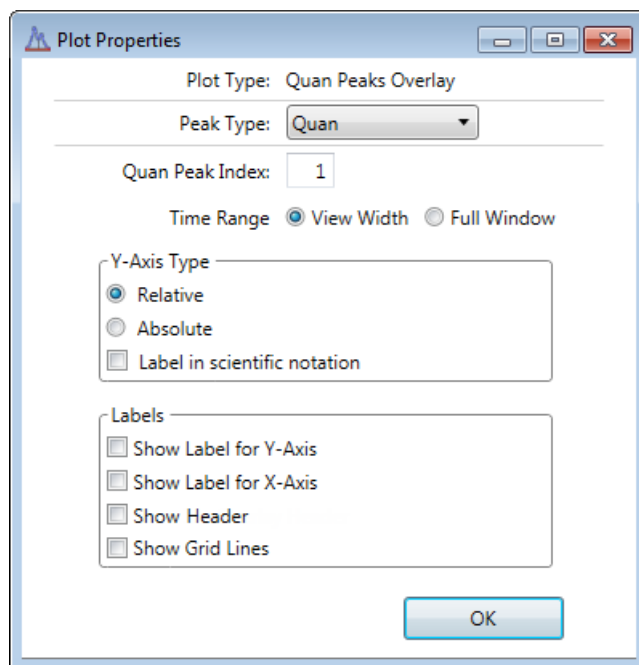


Figure 175. Plot Properties dialog box for a Quan Peaks Overlay plot**Table 89.** Plot Properties parameters for a Quan Peaks Overlay plot

Command	Description
Peak Type	Specifies the type of peak to display in the plot: Quan or InternalStandard
Quan Peak Index	Valid values: 1 through 9
Time Range	Specifies that the plot display one of the following: <ul style="list-style-type: none"> View Width: Zooms the display on the peak of interest. Full Window: Displays the full width of the scan.
Y-Axis Type	
Relative	Displays the <i>y</i> -axis scale from 0 through 100.
Absolute	Displays the <i>y</i> -axis scale from 0 to the actual value of the most intense peak in the group.
Label in Scientific Notation	Displays the <i>y</i> -axis scale in scientific notation.
Labels	
Show Label for Y-Axis	Displays the relative intensity value on the <i>y</i> axis.
Show Label for X-Axis	Displays the retention time (RT) value on the <i>x</i> axis.
Show Header	Displays the raw data file name and compound name.
Show Grid Lines	Displays the grid lines in the plot background.

Spectral Plot

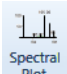
Use the  icon to insert a Spectral plot.

Figure 176. Spectral plot

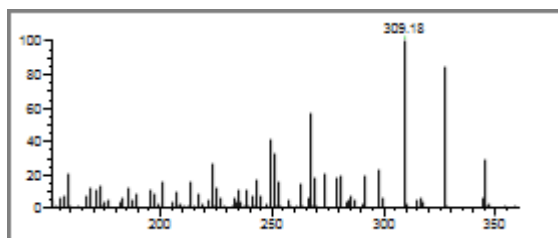


Figure 177. Plot Properties dialog box for a Spectral plot

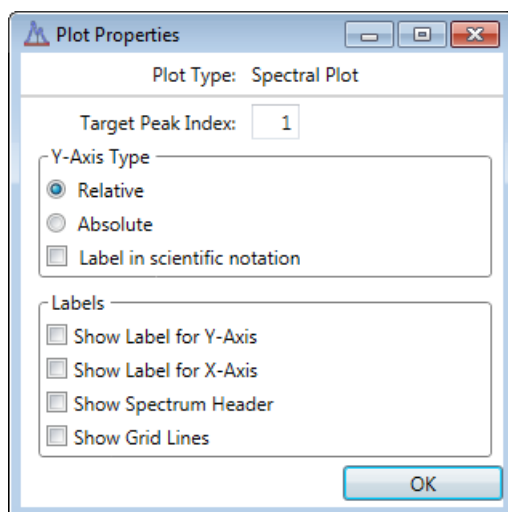


Table 90. Plot Properties parameters for a Spectral plot

Command	Description
Target Peak Index	Valid values: 1 through 9
Y-Axis Type	
Relative	Displays the <i>y</i> -axis scale from 0 through 100.
Absolute	Displays the <i>y</i> -axis scale from 0 to the actual value of the most intense peak in the group.
Label in Scientific Notation	Displays the <i>y</i> -axis scale in scientific notation.
Labels	
Show Label for Y-Axis	Displays the relative intensity value on the <i>y</i> axis.
Show Label for X-Axis	Displays the retention time (RT) value on the <i>x</i> axis.
Show Header	Displays the raw data file name and compound name.
Show Grid Lines	Displays the grid lines in the plot background.

Comparative Spectral Plot (quantitation reports only)


Use the  icon to insert a Comparative Spectral plot.

Figure 178. Comparative Spectral plot

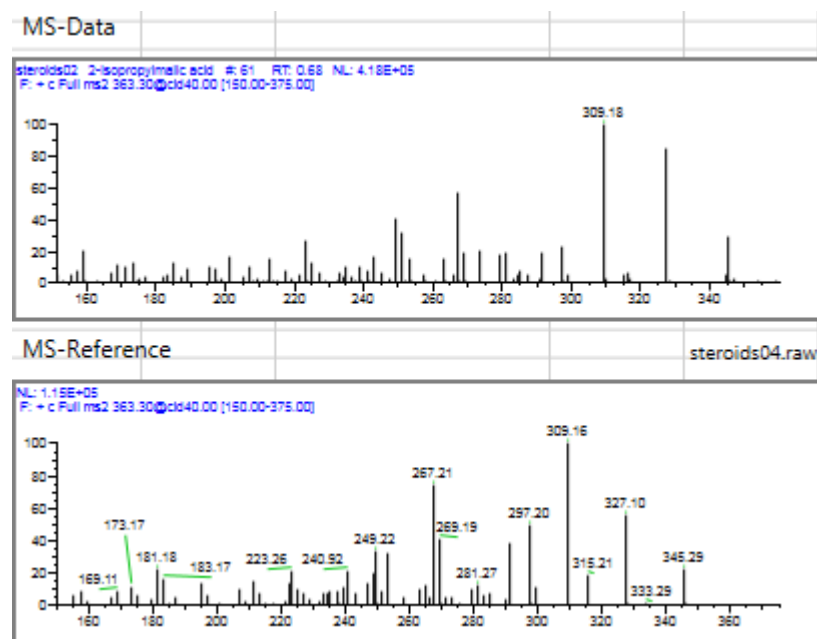


Figure 179. Plot Properties dialog box for a Comparative Spectral plot

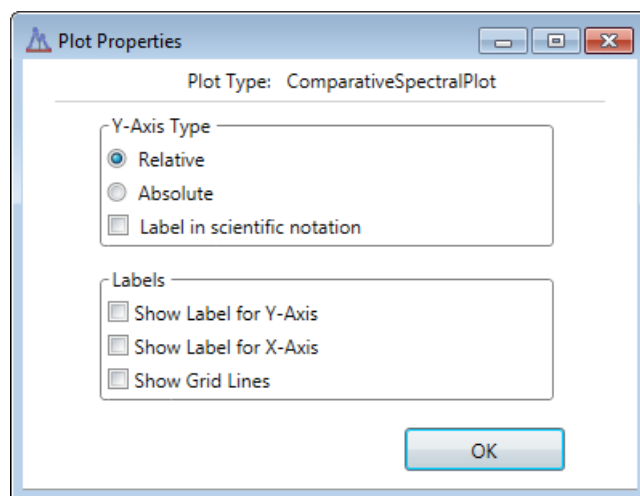


Table 91. Plot Properties parameters for a Comparative Spectral plot (Sheet 1 of 2)

Command	Description
Y-Axis Type	
Relative	Displays the y-axis scale from 0 through 100.

Table 91. Plot Properties parameters for a Comparative Spectral plot (Sheet 2 of 2)

Command	Description
Absolute	Displays the <i>y</i> -axis scale from 0 to the actual value of the most intense peak in the group.
Label in Scientific Notation	Displays the <i>y</i> -axis scale in scientific notation.
Labels	
Show Label for Y-Axis	Displays the relative intensity value on the <i>y</i> axis.
Show Label for X-Axis	Displays the retention time (RT) value on the <i>x</i> axis.
Show Grid Lines	Displays the grid lines in the plot background.

Calibration Curve Plot (quantitation reports only)

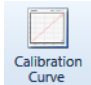
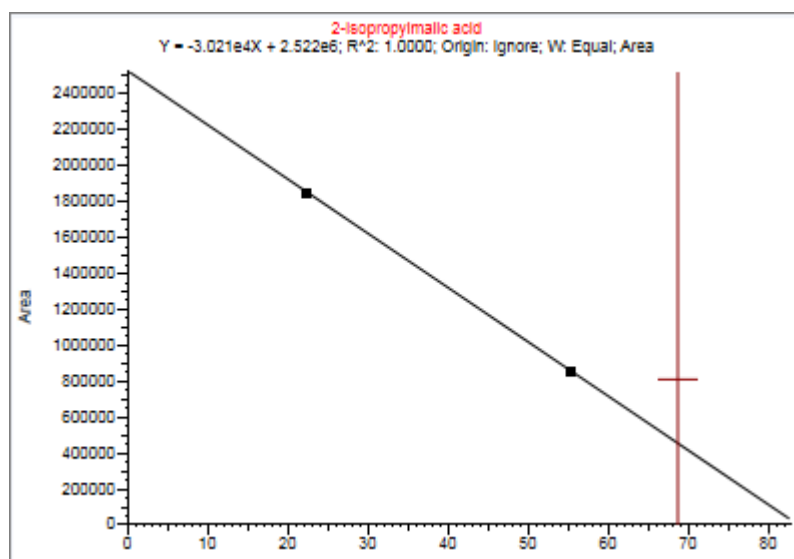
Use the  icon to insert a Calibration Curve plot.


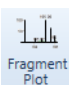

Figure 180. Calibration Curve plot



Note There are no plot properties for a calibration curve plot.

Isotope Plot, Fragment Plot, or Library Plot

Do one of the following:

- Use the  icon to insert an Isotope plot.
- Use the  icon to insert a Fragment plot.
- Use the  icon to insert a Library plot.

The Isotope, Fragment, and Library plots use identical plot properties.

Figure 181. Isotope plot

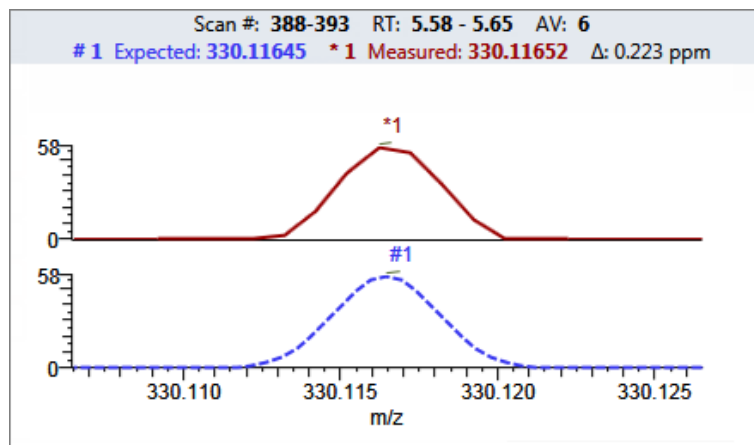


Figure 182. Fragment plot

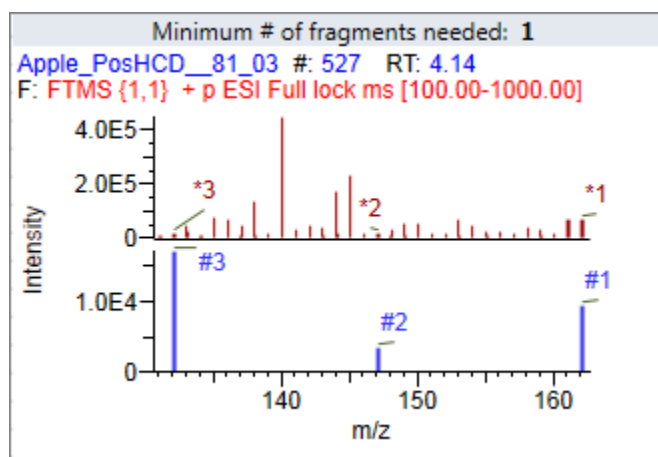


Figure 183. Library plot

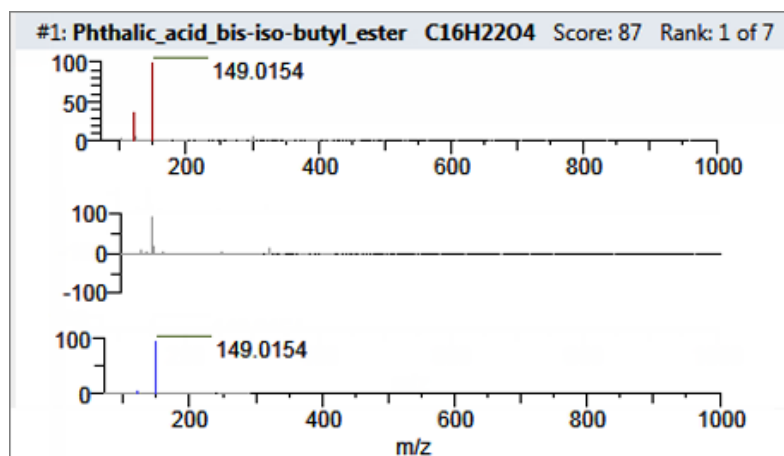


Figure 184. Plot Properties dialog box for an Isotope plot, a Fragment plot, or a Library plot

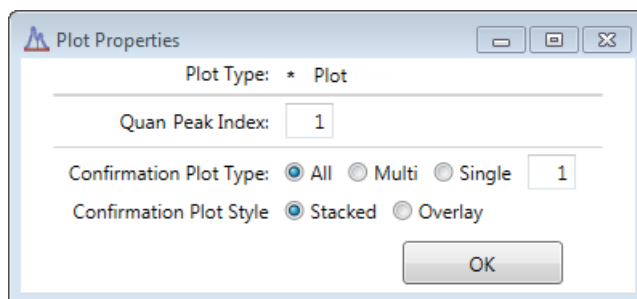


Table 92. Plot Properties parameters for an Isotope plot, a Fragment plot, or a Library plot (Sheet 1 of 2)

Command	Description
Target Peak Index	Specifies the number of peak plots to display. Valid values: 1 through 9
Confirmation Plot Type	Specifies that the plot display one of the following: <ul style="list-style-type: none"> All: Displays isotopes, fragments, or library hits. Multi: Displays individual plots for each isotope. You can individually stack or overlay the plots for each isotope. Single: Displays only the selected isotope, fragment, or library hit. <p>For examples of these display types, see Isotope, Fragments, or Library Match.</p>

Table 92. Plot Properties parameters for an Isotope plot, a Fragment plot, or a Library plot
(Sheet 2 of 2)

Command	Description
Confirmation Plot Style	Specifies that the plot display one of the following: <ul style="list-style-type: none"> Stacked: Stacks the simulated spectrum and the peak apex spectrum. Overlay: Overlays the two spectrum displays. <p>For examples of these display types, see Isotope, Fragments, or Library Match.</p>

Group Bar Graph Plot


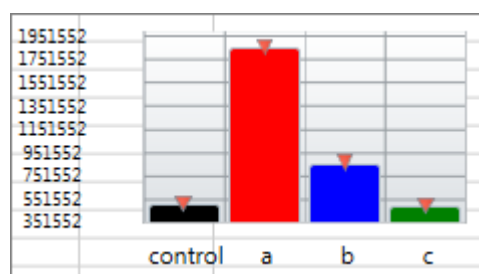
Use the  icon to insert a Group Bar Graph plot.

Figure 185. Group Bar Graph plot



Note There are no plot properties for a group bar graph plot.

RT Summary Plot


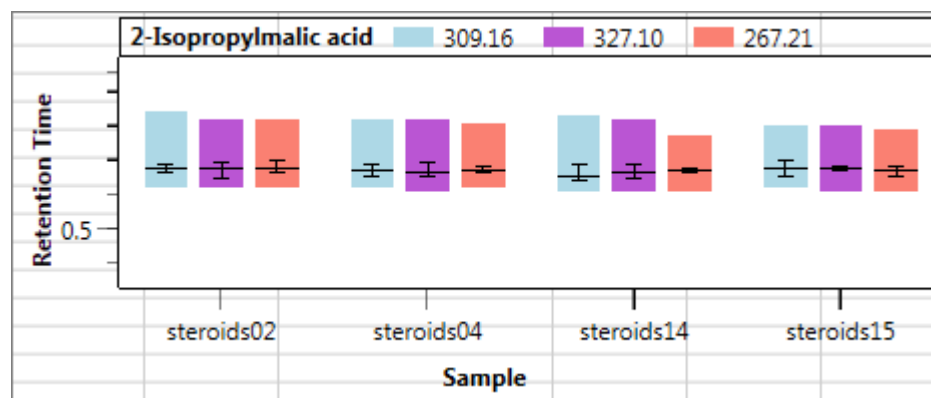
Use the  icon to insert an RT Summary plot.

Figure 186. RT Summary plot



Note There are no plot properties for an RT summary plot.

XIC Overlay Plot

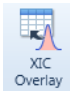
Use the  icon to insert an XIC Overlay plot.

Figure 187. XIC Overlay plot

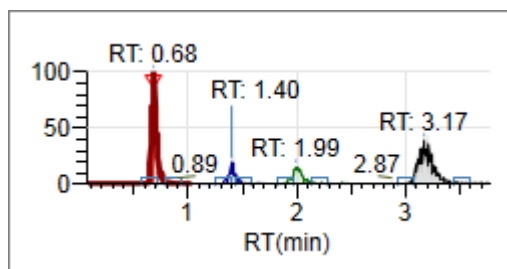


Figure 188. Plot Properties dialog box for an XIC Overlay plot

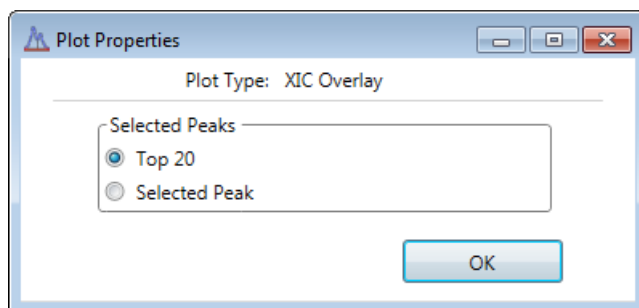


Table 93. Plot Properties parameters for an XIC Overlay plot

Command	Description
Selected Peaks	
Top 20	Displays the top 20 peaks.
Selected Peak	Displays only the selected peak.

Batch Quan Peaks Overlay Plot

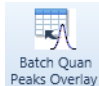
Use the  icon to insert a Batch Quan Peaks Overlay plot.

Figure 189. Batch Quan Peaks Overlay plot

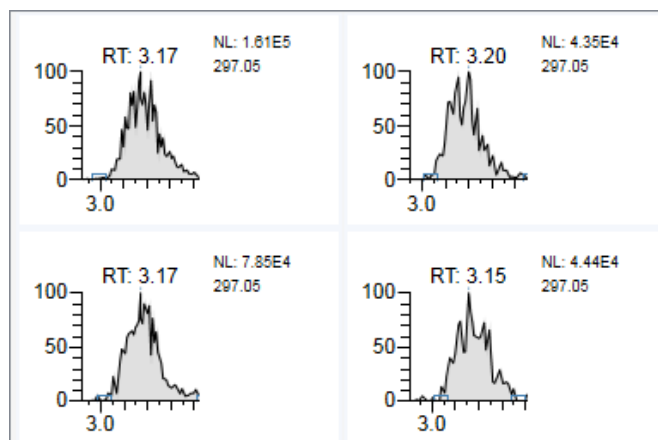


Figure 190. Plot Properties dialog box for a Batch Quan Peaks Overlay plot

Table 94. Plot Properties parameters for a Batch Quan Peaks Overlay plot (Sheet 1 of 2)

Command	Description
Quan Peak Index	Valid values: 1 through 9
Time Range	Specifies that the plot display one of the following: <ul style="list-style-type: none"> View Width: Zooms the display on the peak of interest. Full Window: Displays the full width of the scan.

Table 94. Plot Properties parameters for a Batch Quan Peaks Overlay plot (Sheet 2 of 2)

Command	Description
Y-Axis Type	
Relative	Displays the <i>y</i> -axis scale from 0 through 100.
Absolute	Displays the <i>y</i> -axis scale from 0 to the actual value of the most intense peak in the group.
Label in Scientific Notation	Displays the <i>y</i> -axis scale in scientific notation.
Labels	
Show Label for Y-Axis	Displays the relative intensity value on the <i>y</i> axis.
Show Label for X-Axis	Displays the retention time (RT) value on the <i>x</i> axis.
Show Header	Displays the raw data file name and compound name.
Show Grid Lines	Displays the grid lines in the plot background.
Number of Columns	Specifies the number of columns that you want to display in the inserted graphic. The application resizes the images to fit the specified number of columns.

Unknown Screening Plots

To customize the display in the unknown screening report, use the Plot Properties dialog box for each plot.

- [Peak Chromatogram Plot](#)
- [TIC Chromatogram Plot](#)
- [Spectral Plot](#)
- [XIC Overlay Plot](#)
- [Library Plot, Fragments Plot, or Isotopes Plot](#)
- [XIC Plot](#)
- [Cross Sample Peak Overlay Plot](#)
- [Chemical Structure Plot](#)
- [Group Bar Graph Plot](#)

Peak Chromatogram Plot

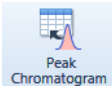
Use the  icon to insert a Peak Chromatogram plot.

Figure 191. Peak Chromatogram plot

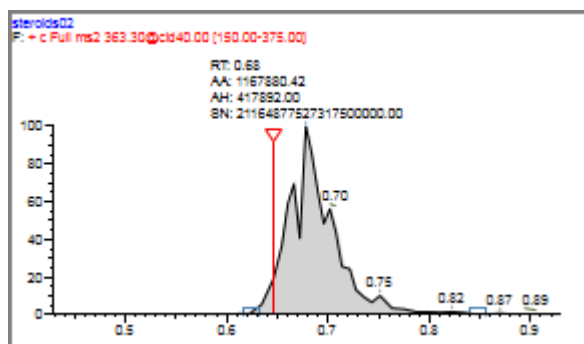


Figure 192. Plot Properties dialog box for Peak Chromatogram plot

Table 95. Plot Properties parameters for a Peak Chromatogram plot (Sheet 1 of 2)

Command	Description
Y-Axis Type	
Relative	Displays the <i>y</i> -axis scale from 0 through 100.
Absolute	Displays the <i>y</i> -axis scale from 0 to the actual value of the most intense peak in the group.
Label in Scientific Notation	Displays the <i>y</i> -axis scale in scientific notation.
Labels	
Show Label for Y-Axis	Displays the relative intensity value on the <i>y</i> axis.

Table 95. Plot Properties parameters for a Peak Chromatogram plot (Sheet 2 of 2)

Command	Description
Show Label for X-Axis	Displays the retention time (RT) value on the <i>x</i> axis.
Show Grid Lines	Displays the grid lines in the plot background.
Peak Plot Labels	
RT	Displays the retention time (RT) label in the peak chromatogram plot.
Area	Displays the peak area (AA) label in the peak chromatogram plot.
Height	Displays the peak height (AH) label in the peak chromatogram plot.
Signal to Noise	Displays the signal-to-noise (SN) label in the peak chromatogram plot.
M/Z	Displays the peak mass.
Show All Apex Times	Displays all surrounding apex times in the plot.
Sample Filename	Displays the name of the sample file.
Filter	Displays the filter used for the scan. The plot displays the filter name below the raw data file name.

TIC Chromatogram Plot

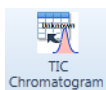
Use the  icon to insert a TIC Chromatogram plot.

Figure 193. TIC Chromatogram plot

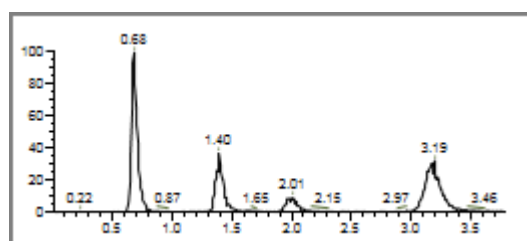


Figure 194. Plot Properties dialog box for a TIC Chromatogram plot

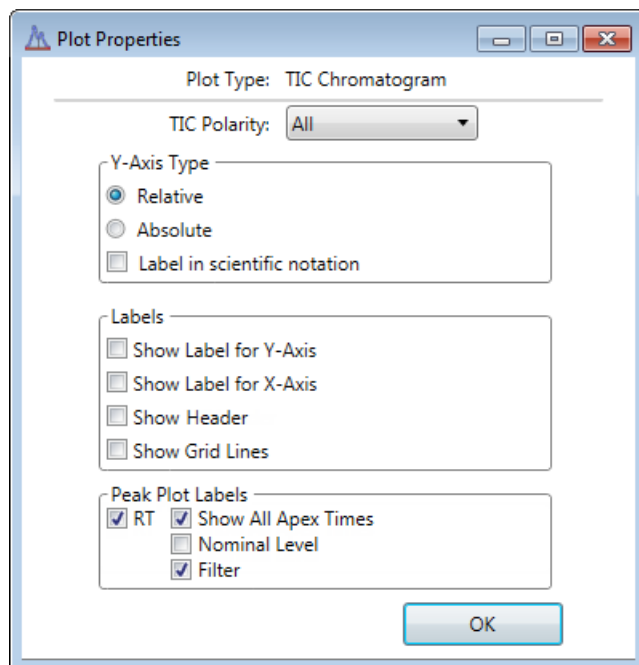


Table 96. Plot Properties parameters for a TIC Chromatogram plot (Sheet 1 of 2)

Command	Description
TIC Polarity	Specifies the polarity of adducts displayed in the plot. All: Displays positive, negative, and analog adducts. Positive: Displays only positive adducts. Negative: Displays only negative adducts. Analog: Displays only analog adducts.
Y-Axis Type	
Relative	Displays the <i>y</i> -axis scale from 0 through 100.
Absolute	Displays the <i>y</i> -axis scale from 0 to the actual value of the most intense peak in the group.
Label in Scientific Notation	Displays the <i>y</i> -axis scale in scientific notation.
Labels	
Show Label for Y-Axis	Displays the relative intensity value on the <i>y</i> axis.
Show Label for X-Axis	Displays the retention time (RT) value on the <i>x</i> axis.
Show Header	Displays the raw data file name.
Show Grid Lines	Displays the grid lines in the plot background.
Peak Plot Labels	
RT	Displays the retention time (RT) label in the TIC chromatogram plot.
Show All Apex Times	Displays all surrounding apex times in the plot.

Table 96. Plot Properties parameters for a TIC Chromatogram plot (Sheet 2 of 2)

Command	Description
Nominal Level	Displays the nominal level (NL) values for each peak in the plot.
Filter	Displays the filter used for the scan. The plot displays the filter name below the raw data file name.

Spectral Plot

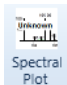
Use the  icon to insert a Spectral plot.

Figure 195. Spectral plot

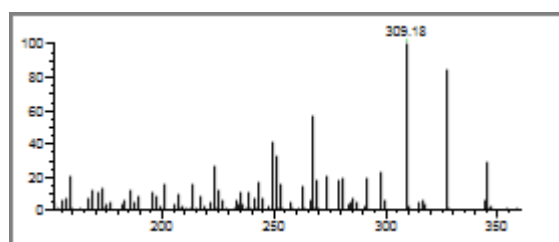


Figure 196. Plot Properties dialog box for a Spectral plot

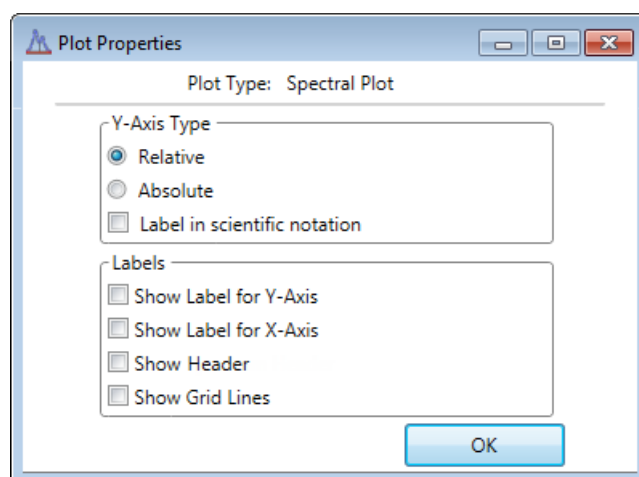


Table 97. Plot Properties parameters for a Spectral plot (Sheet 1 of 2)

Command	Description
Y-Axis Type	
Relative	Displays the y-axis scale from 0 through 100.
Absolute	Displays the y-axis scale from 0 to the actual value of the most intense peak in the group.
Label in Scientific Notation	Displays the y-axis scale in scientific notation.

Table 97. Plot Properties parameters for a Spectral plot (Sheet 2 of 2)

Command	Description
Labels	
Show Label for Y-Axis	Displays the relative intensity value on the <i>y</i> axis.
Show Label for X-Axis	Displays the retention time (RT) value on the <i>x</i> axis.
Show Header	Displays the name of the sample file, the name of the compound, the retention time, the nominal level, and the filter used for the scan.
Show Grid Lines	Displays the grid lines in the plot background.

XIC Overlay Plot

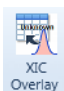
Use the  icon to insert an XIC Overlay plot.

Figure 197. XIC Overlay plot

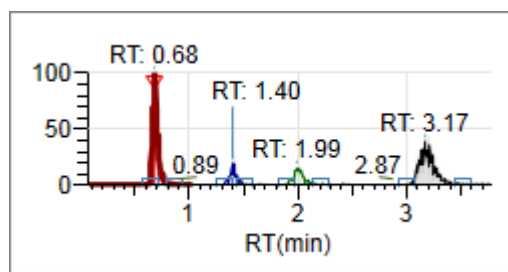


Figure 198. Plot Properties dialog box for an XIC Overlay plot

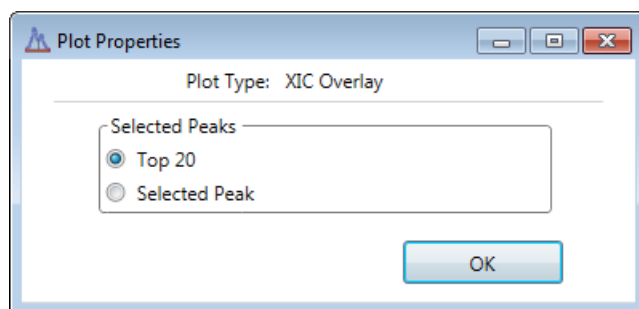
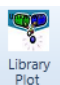
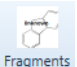



Table 98. Plot Properties parameters for an XIC Overlay plot

Command	Description
Selected Peaks	
Top 20	Displays the top 20 peaks.
Selected Peak	Displays only the selected peak.

Library Plot, Fragments Plot, or Isotopes Plot

Do one of the following:

- Use the  icon to insert a Library plot.
- Use the  icon to insert a Fragments plot.

Use the  icon to insert an Isotope plot.

The Library, Fragments, and Isotopes plots use identical plot properties.

Figure 199. Library plot

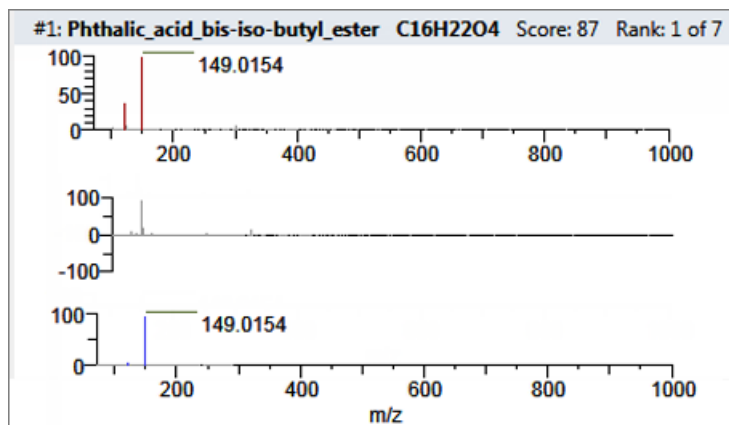


Figure 200. Fragments plot

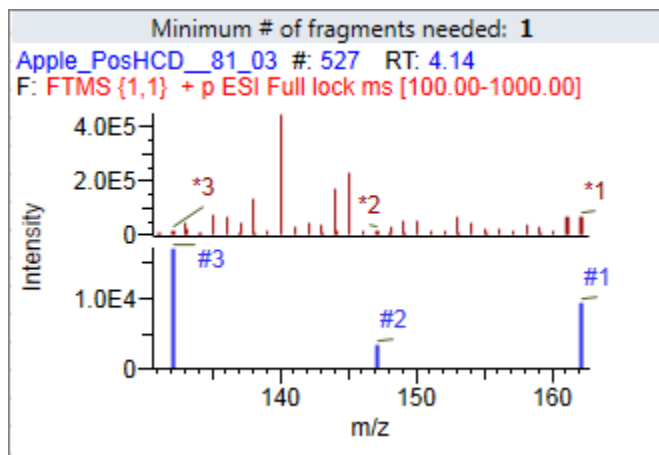


Figure 201. Isotopes plot

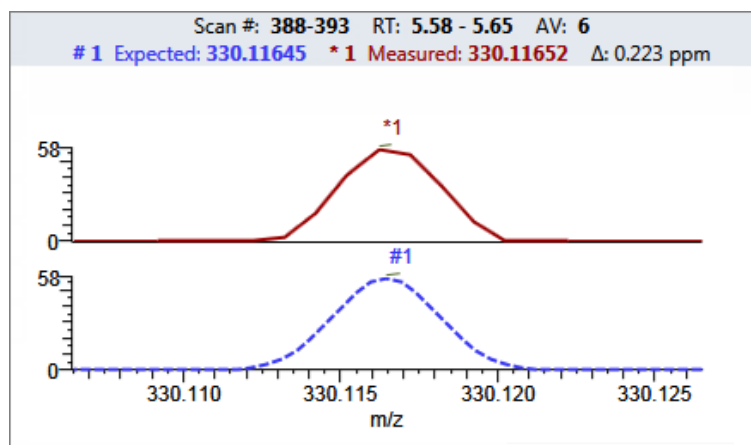


Figure 202. Plot Properties dialog box for a Library plot, a Fragments plot, or an Isotopes plot

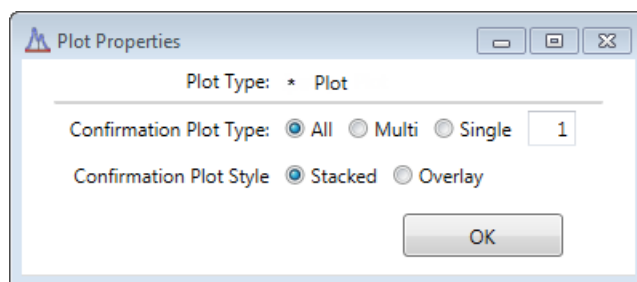


Table 99. Plot Properties parameters for a Library plot, a Fragments plot, or an Isotopes plot

Command	Description
Confirmation Plot Type	<p>Specifies that the plot display one of the following:</p> <ul style="list-style-type: none"> All: Displays isotopes, fragments, or library hits. Multi: Displays individual plots for each isotope. You can individually stack or overlay the plots for each isotope. Single: Displays only the selected isotope, fragment, or library hit. <p>For examples of these display types in the Data Review, see Library Search Pane, Fragments Pane, or Isotopes Pane.</p>
Confirmation Plot Style	<p>Specifies that the plot display one of the following:</p> <ul style="list-style-type: none"> Stacked: Stacks the simulated spectrum and the peak apex spectrum. Overlay: Overlays the two spectrum displays. <p>For examples of these display types in the Data Review, see Library Search Pane, Fragments Pane, or Isotopes Pane.</p>

XIC Plot

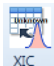
Use the  icon to insert an XIC plot.

Figure 203. XIC plot

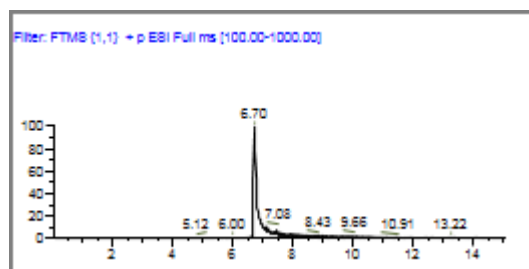


Figure 204. Plot Properties dialog box for an XIC plot

Table 100. Plot Properties parameters for an XIC plot (Sheet 1 of 2)

Command	Description
TIC Polarity	Specifies the polarity of adducts displayed in the plot. All: Displays positive, negative, and analog adducts. Positive: Displays only positive adducts. Negative: Displays only negative adducts. Analog: Displays only analog adducts.
Y-Axis Type	
Relative	Displays the y-axis scale from 0 through 100.

Table 100. Plot Properties parameters for an XIC plot (Sheet 2 of 2)

Command	Description
Absolute	Displays the y -axis scale from 0 to the actual value of the most intense peak in the group.
Label in Scientific Notation	Displays the y -axis scale in scientific notation.
Labels	
Show Label for Y-Axis	Displays the relative intensity value on the y axis.
Show Label for X-Axis	Displays the retention time (RT) value on the x axis.
Show Header	
Show Grid Lines	Displays the grid lines in the plot background.
Peak Plot Labels	
RT	Displays the retention time (RT) values for each peak in the plot.
Show All Apex Times	Displays all surrounding apex times in the plot.
Nominal Level	Displays the nominal level (NL) values for each peak in the plot.
Filter	Displays the filter used for the scan. The plot displays the filter name below the raw data file name.

Cross Sample Peak Overlay Plot

Use the  icon to insert a Cross Sample Peak Overlay plot.

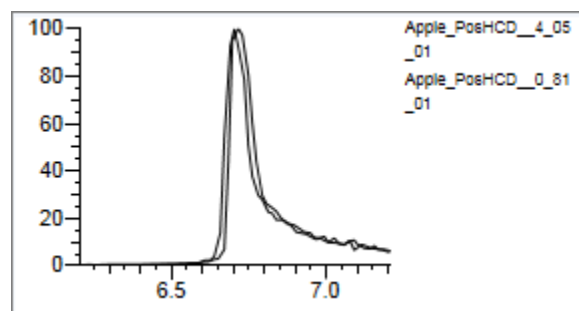
Figure 205. Cross Sample Peak Overlay plot


Figure 206. Plot Properties dialog box for a Cross Sample Peak Overlay plot

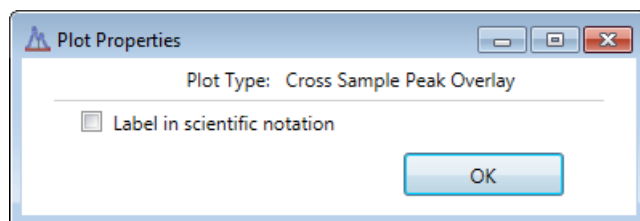


Table 101. Plot Properties parameter for a Cross Sample Peak Overlay plot

Command	Description
Label in Scientific Notation	Displays the y-axis scale in scientific notation.

Chemical Structure Plot

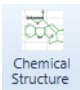
Use the  icon to insert a Chemical Structure plot.

Figure 207. Chemical Structure plot

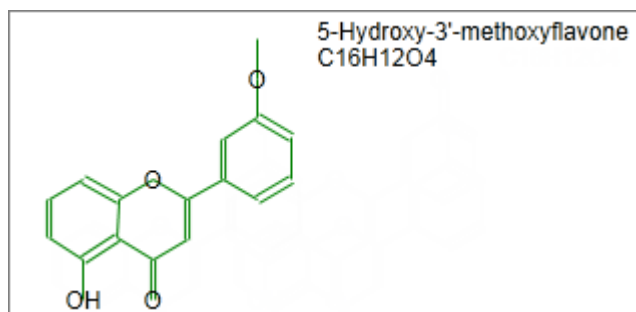


Figure 208. Plot Properties dialog box for a Chemical Structure plot

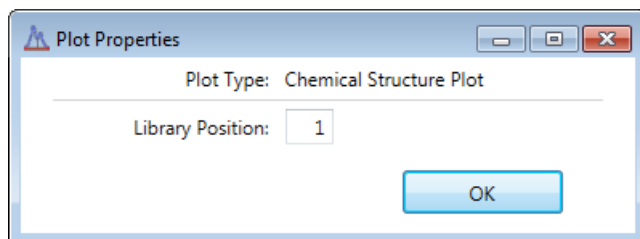


Table 102. Plot Properties parameter for a chemical structure plot

Command	Description
Library Position	Specifies the hierarchical position in the library when there are multiple chemical structures.

Group Bar Graph Plot


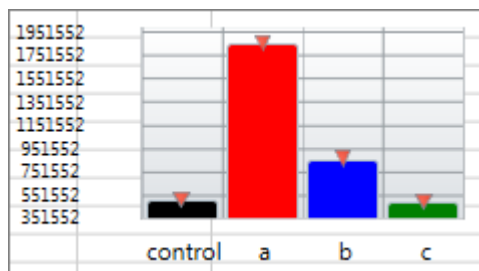
Use the  icon to insert a Group Bar Graph plot.

Figure 209. Group Bar Graph plot



Note There are no plot properties for a group bar graph plot.

Quick Tips

Follow these tips to make the most of the Report Designer features.

❖ To move a report to the Archive folder

1. With a report open for editing, choose **File > Archive** from the Report Designer main menu.
2. At the confirmation prompt, click **Yes**.

The application deletes the report from the ...*batch*\ReportOutput folder.

❖ To insert items in the Report Designer

- Insert a data table when the selected cell is above or below any other table, but not in the same row as a field. You cannot insert tables in repeat areas.
- Insert a data field when the selected cell is not in the same row as a table.
- Use these shortcut keys:

CTRL+T	Insert a table.
CTRL+T	Edit a table when the selected cell is inside a table.
CTRL+SHIFT+T	Insert or edit a field.
CTRL+R	Insert or edit a repeating area.

❖ To format cells and group headers

Because the first row of a data table retains formatting information, edit the formatting of the first row.

The application copies the formatting to all other rows.

For group header items, the application copies the formatting of the first group header to the remaining group headers.

❖ To sort fields

Select a field, and then click **A-Z** to sort the data in ascending or descending order.

Using the Audit Viewer

The TraceFinder application records all user access, including logging in, logging out, data creation and editing (batches, methods, and templates), and manual integration. You can use the [Audit Viewer](#) to view the resulting log files to track modifications to the data. When an event requires confirmation (as specified in the Administrator Console), the Audit Viewer records who confirmed each change to a batch, method, or template. When no confirmation is required, then the Audit Viewer records the user who was logged in when the change occurred.

IMPORTANT TraceFinder 4.1 uses the same data as TraceFinder 4.0. By default, the application stores the log files for the 4.1 release in the TraceFinderData\4.0 or TraceFinder\4.0 folders.

In the Administration Console, a user with Auditing permissions can configure the auditing service by specifying which events are logged, which events require confirmation, a list of default reasons for a specific event, and whether a user can submit a custom reason. To use the auditing administration tools, refer to the instructions in the *TraceFinder Administrator Console User Guide*.

Contents

- [Audit Trail Log Files](#)
- [Audit Log File Functions](#)
- [Audit Viewer](#)

Audit Trail Log Files

The application creates the following audit trail log files:

- **Application:** Records all user access, such as starting and stopping the application, logging in, logging out, or accessing or saving data in batches and methods. The application saves the data in the following log file: C:\Thermo\TraceFinder\4.0\Logs\AuditLog.adb.
- **Master Method:** Records all user interactions with master methods, such as creating, opening, or editing a master method. The application saves the data in the following log file: C:\TraceFinderData\4.0\Methods*MasterMethodName*\AuditLog.adb.

- **Batch Template:** Records all user interactions with batch templates, such as creating, opening, or editing a batch template. The application saves the data in the following log file: C:\TraceFinderData\4.0\Templates\Batches\BatchTemplateName\AuditLog.adb.
- **Batch:** Records all user interactions with batches, such as creating, opening, editing, acquiring, processing, or generating reports for a batch. The application saves the data in the following log file:
C:\TraceFinderData\4.0\Projects\SubFolder\BatchName\AuditLog.adb.

Audit Log File Functions


The Audit Viewer displays all saved audit log files, and you can filter and sort the audit data.

Note The Audit Viewer is available only when you have enabled auditing from the Administration Console. Refer to the *TraceFinder Administrator Console User Guide*.

Follow these procedures:

- [To access the Audit Viewer](#)
- [To select an audit log](#)
- [To view only application, method, batch, or batch template events](#)
- [To create a filter for audit log events](#)
- [To view audit event details](#)
- [To create a filter for an audit log history](#)
- [To display the history for an event](#)

❖ To access the Audit Viewer

Choose **Tools > Audit Trail** from the TraceFinder main menu or click the **Audit Viewer** icon,  .

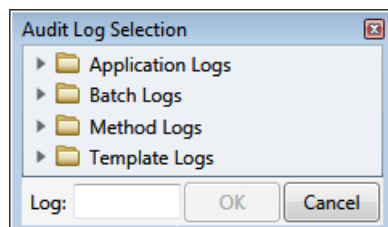
The [Audit Viewer](#) opens.

Note The Tools > Audit Trail menu command always opens to application log files, whereas the Audit Viewer icon is context sensitive and opens to the appropriate type of log files (application, method, batch, or batch template).

❖ **To select an audit log**

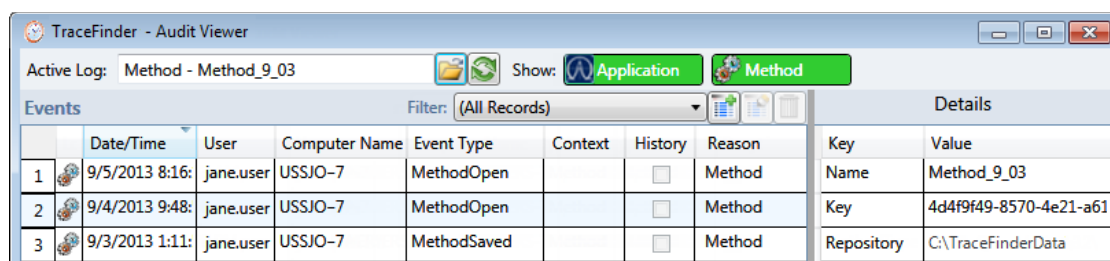
1. Click the **Open Audit Log** icon, .

The application opens the Audit Log Selection dialog box.



2. Expand a log folder to select an application, batch, method, or batch template audit log file, and click **OK**.

The Audit Viewer displays the contents of the selected audit log file, as in this example for a method.



❖ **To view only application, method, batch, or batch template events**

1. In the Active Log list, select an Active Log file.
 - Application logs include application and security events.
 - Method logs include application and method events.
 - Batch logs include application, method, and batch events.
 - Batch template logs include application and batch template events.

The Audit Viewer displays icons for each of the event types included in the log file.



2. Click an icon to turn the display on or off.

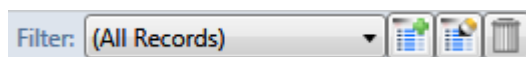
In this example, the Events pane displays Batch and Method events and hides Application events.



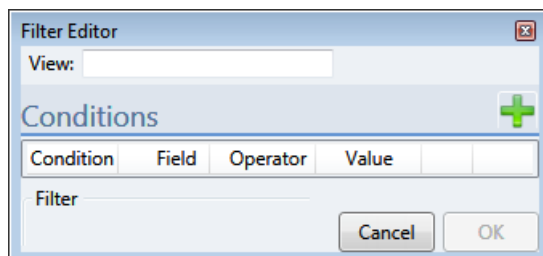
Note There is no icon for security events in an application log file. You cannot hide the display of security events.

❖ **To create a filter for audit log events**

1. In the Events pane, click the **Create New Filter** icon, .



The Filter Editor dialog box opens.

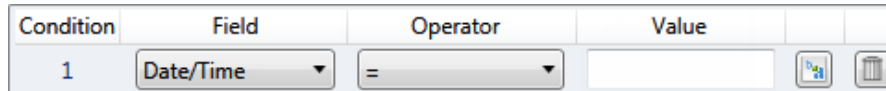


2. In the View box, type a name for the new filter.

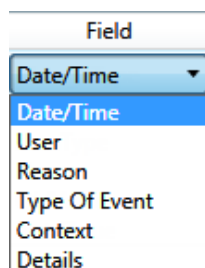
You can also leave the View box empty when you enter your filter criteria. When you finish adding conditions and click OK (step 8), the application filters the current events list based on the filter criteria you specify, but the filter is not saved. The Filter list in the viewer identifies this filter as (Custom).

3. Click the **Add** icon, .

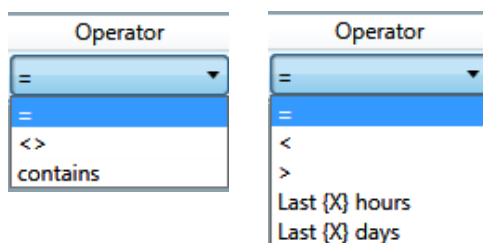
The application adds a new, undefined condition to the Condition list.



4. In the Field list, select one of the following field types.

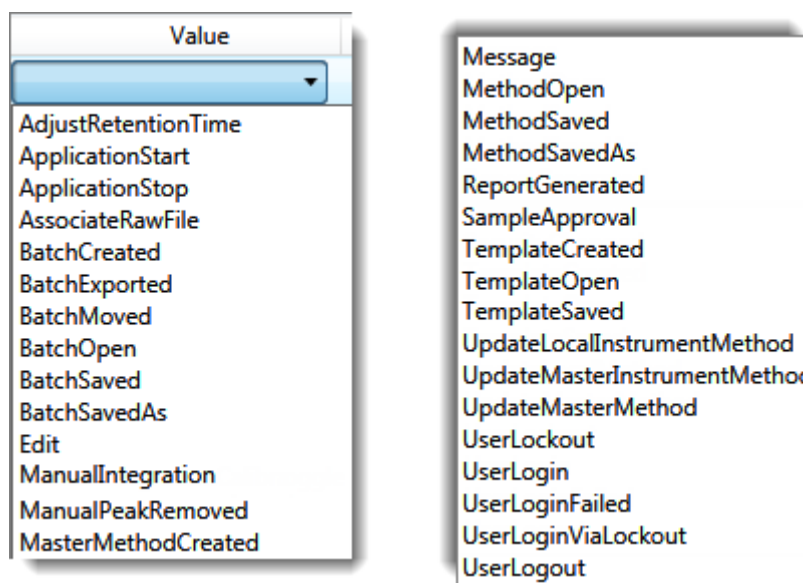



5. In the Operator list, select one of the available operators.



The available operators depend on the field type that you selected.

6. In the Value box, type a value or select a value from the list.
 - For the Date/Time field, type a numerical value in the Value box.
 - For the User, Reason, Context, or Details field, type the appropriate text in the Value box. This value is case sensitive.
 - For the Type of Event field, select one of the following values.



Note You can also click the Tokens icon, , and select a token for the Value. Tokens are predefined values, such as dates and sample or compound identifiers.

Condition	Field	Operator	Value
1	Date/Time	=	{Today}

Filter
((TimeStamp >= datetime("2013-11-18T00:00:00") and TimeStamp <= datetime("2015-01-18T23:59:59")))

The Filter at the bottom of the dialog box displays the complete definition for the filter.

Condition	Field	Operator	Value
1	User	=	jane.user

Filter
(((Userid like '%\jane.user') or (Userid = 'jane.user')))

7. Repeat [step 3](#) (on [page 434](#)) through [step 6](#) (on this page) for each condition that you want to include in your filter.
8. When you have added all your conditions, click **OK**.

The application creates the new filter with the specified conditions.

❖ To view audit event details

In the Events pane, select an event.

The Details pane displays key values based on the type of log you select.



Details for batch log files include the name of the batch.



Details for method log files include the name of the method and the method type.



Details for batch template log files include the name of the batch template, the location of the data repository, and the subproject folder where the template was created.




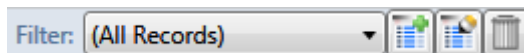
Details for Application log files include whether the events are for a batch or method, the location of the data repository, and, for a batch, the subproject folder where the batch was created.



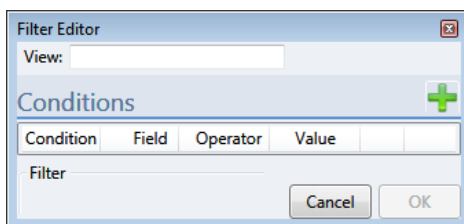
Details for Security log files include the authentication method used (Windows Active Directory or local machine) and the location of the administrator repository.

❖ To create a filter for an audit log history

1. In the History pane, click the **Create New Filter** icon, .



The Filter Editor dialog box opens.

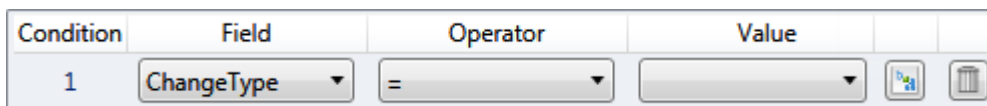


2. In the View box, type a name for the new filter.

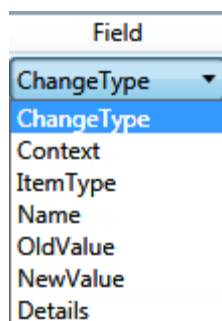
You can also leave the View box empty when you enter your filter criteria. When you finish adding conditions and click OK (step 8), the application filters the current history list based on the criteria you specify, but the filter is not saved. The Filter list in the viewer identifies this filter as (Custom).

3. Click the **Add** icon, .

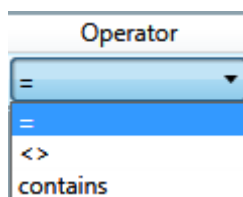
The application adds a new, undefined condition to the Conditions list.



4. In the Field list, select one of the following field types.

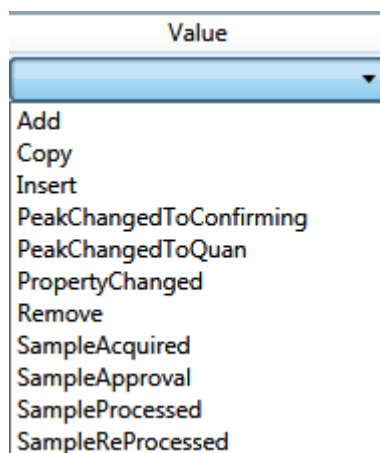


5. In the Operator list, select from *equals*, *less than/greater than*, or *contains*.



The available operators depend on the field type that you selected.

6. In the Value list, type or select a value from the list.
- For the Name, Context, or Details field, type the appropriate text in the Value box. This value is case sensitive.
 - For the OldValue or NewValue field, type a numerical value in the Value box.
 - For the ChangeType field, select one of the following values.



- When the selected Field is **ItemType**, select one of the following values.

Value

Application

AutoSample

Batch

BatchTemplate

BreakDownGroup

CalibrationHistory

CalibrationLevel

Compound

CompoundCalLevel

CompoundGroup

CompoundQCLLevel

CompoundTransition

DetectionParms

InstrumentMethod

IntelligentSequence

MassRange

Method

PeakIdentifier

QCLLevel

ReferenceSample

Role

RoleMap

RolePermission


Sample



Security

Suitability

User

UserRole



Note You can also click the Tokens icon, , and select a token for the Value. Tokens are predefined values, such as dates and sample or compound identifiers.

Condition	Field	Operator	Value		
1	Name	=	{User}		

Filter

((Name = 'jane.user'))

The Filter at the bottom of the dialog box displays the complete definition for the filter.

Condition	Field	Operator	Value		
1	ChangeType	=	PropertyChanged		

Filter

((ChangeType = 'PropertyChanged'))

- Repeat steps 3 through 6 for each condition that you want to include in your filter.
- When you have added all your conditions, click **OK**.

The applications created the new filter with the specified conditions.

❖ To display the history for an event

- In the Events pane, select an event that has a selected check box in the History column.

Event Type	Context	History	Reason
BatchOpen		<input type="checkbox"/>	Batch
BatchCreated		<input type="checkbox"/>	Batch
BatchSaved		<input checked="" type="checkbox"/>	Batch

Select an event that has the History check box selected.

The History pane displays all unsaved (queued) actions for the selected event.

History							Filter: (All Records)
	Order	Change Type	Context	Item Type	Item Name	Old Value	New Value
1	1	Add	Unknown2	Sample	n/a	n/a	1
2	2	Add	Unknown3	Sample	n/a	n/a	2
3	3	Add	Unknown4	Sample	n/a	n/a	3
4	4	PropertyChanged	Apple_0_81	Sample	RawFileName	Unknown1	Apple_0_81_01
5	5	PropertyChanged	Apple_0_81	Sample	SampleID		1
6	6	PropertyChanged	Apple_0_81	Sample	VialPosition		CStk1-01:14
7	7	PropertyChanged	Apple_0_81	Sample	BarcodeExpected		2134

- To limit the actions in the history list, select a filter from the Filter list.

See [To create a filter for an audit log history](#).

Figure 210. Example History filter

When you select this filter,

The Filter Editor dialog box is shown with the following details:

- View:** Column Value Changes
- Conditions:** A table with one condition:

Condition	Field	Operator	Value
1	ChangeType	=	PropertyChanged
- Filter:** (Empty text field)
- Buttons:** OK, Cancel

the History pane displays only these actions for the selected event:

History							Filter: Column Value Changes
	Order	Change Type	Context	Item Type	Item Name	Old Value	New Value
1	1	PropertyChanged	Apple_0_81	Sample	RawFileName	Unknown1	Apple_0_81_01
2	2	PropertyChanged	Apple_0_81	Sample	SampleID		1
3	3	PropertyChanged	Apple_0_81	Sample	VialPosition		CStk1-01:14
4	4	PropertyChanged	Apple_0_81	Sample	BarcodeExpected		2134

Audit Viewer

Use the Audit Viewer to view the audit log files to track user access and modifications to the data. See [Audit Viewer parameters](#).

Figure 211. Application log in the Audit Viewer

TraceFinder - Audit Viewer

Active Log: Application - Logs Show: Application

Events Filter: (All Records)

Log	Date/Time	User	Computer Name	Event Type	Context	History	Reason
1	10/24/2013 1:2	jane.user	USSJO-7	UserLogin		<input type="checkbox"/>	Security
2	10/24/2013 1:2		USSJO-7	ApplicationStart	1.0	<input type="checkbox"/>	Application
3	10/24/2013 11	jane.user	USSJO-7	ApplicationStop		<input type="checkbox"/>	Security
4	10/24/2013 11		USSJO-7	ApplicationStart		<input type="checkbox"/>	Security
5	10/24/2013 10	jane.user	USSJO-7	ApplicationStop		<input type="checkbox"/>	Application
6	10/24/2013 10	jane.user	USSJO-7	UserLogin		<input type="checkbox"/>	Security
7	10/24/2013 10		USSJO-7	ApplicationStart	1.0	<input type="checkbox"/>	Application

Details

Key	Value
UserId	jane.user
UserIsValid	True
UserIsEnabled	True
SecurityEnabled	Yes
SecurityMode	ActiveDirectory
SecurityDomain	AMER
FromLockout	False

History Filter: (All Records)

Order	Change Type	Context	Item Type	Item Name	Old Value	New Value
-------	-------------	---------	-----------	-----------	-----------	-----------

Details

Key	Value
UniqueId	4d5a1806-4f23-4f0

Logged in user: AMER\jane.user

Figure 212. Batch log in the Audit Viewer

TraceFinder - Audit Viewer

Active Log: Batch - Project1\Apple_10_17 Show: Application Batch Method

Events Filter: (All Records)

Log	Date/Time	User	Computer Name	Event Type	Context	History	Reason
1	10/24/2013 8:53	jane.user	USSJO-7	BatchOpen	Apple_10_17	<input type="checkbox"/>	Batch
2	10/24/2013 7:31	jane.user	USSJO-7	BatchSaved	Apple_10_17	<input checked="" type="checkbox"/>	Batch
3	10/24/2013 6:51	jane.user	USSJO-7	BatchSaved	Apple_10_17	<input checked="" type="checkbox"/>	Batch

Details

Key	Value
Repository	C:\TraceFinderData
SubFolder	Project1

History Filter: (All Records)

Order	Change Type	Context	Item Type	Item Name	Old Value	New Value
1	1	Add	Unknown2	Sample	n/a	1
2	2	Add	Unknown3	Sample	n/a	2
3	3	Add	Unknown4	Sample	n/a	3
4	4	PropertyChanged	Apple_0_81_01	Sample	RawFileName	Unknown1
5	5	PropertyChanged	Apple_0_81_01	Sample	SampleID	1
6	6	PropertyChanged	Apple_0_81_01	Sample	VialPosition	CSstk-01:14

Details

Key	Value
UniqueId	4d5a1806-4f23-4f03

Logged in user: AMER\jane.user

Table 103. Audit Viewer parameters (Sheet 1 of 2)
















Parameter	Description
Active Log	Name of the current audit log file.
	Opens the Audit Log Selection dialog box where you can open a different audit log file. You can select from these audit log files: application, batch, method, or batch template.
	Refreshes the current audit log file in the viewer.
Show	Select to display only specific types of events in the Events list. The selected log file can contain batch, method, and application events.
Events	
Filter	<p>Select a filter view to use for displaying the event log entries.</p>  <p> Opens the Filter Editor dialog box where you can create a filter view.</p> <p> Opens the Filter Editor dialog box where you can edit the current filter view.</p> <p> Deletes the current event filter view.</p>
Log	<p> Indicates an event that occurred at the main application level, such as logging in or opening a batch.</p> <p> Indicates an event that occurred in the Administration Console.</p> <p> Indicates an event that occurred in a method template.</p> <p> Indicates an event that occurred in a method.</p> <p> Indicates an event that occurred in a batch.</p>
Date/Time	Time stamp of the event.
User	<p>When an event requires confirmation (as specified in the Administrator Console), <i>User</i> is the user who confirmed each change to a batch, method, or template.</p> <p>When no confirmation is required, <i>User</i> is the user who was logged in when the change occurred.</p>
Computer Name	Name of the computer on which the application recorded the event.
Event Type	Specific event that triggered the log file entry. For a complete list of event types, refer to the <i>TraceFinder Administration Console User Guide</i> .
Context	Name of the sample, batch, method, or application version where the event occurred.

Table 103. Audit Viewer parameters (Sheet 2 of 2)


Parameter	Description
History	Indicates that there is a history log of queued actions for the event.
Reason	Default or custom reason that the user entered for the event.
Details Key/Value	Identifying parameters and their values for the selected auditing event. These key parameters are different for each type of auditing event.
History	
Filter	<p>Select a filter view to use for displaying the change history.</p>  <p>  Opens the Filter Editor dialog box where you can create a filter view. </p> <p>  Opens the Filter Editor dialog box where you can edit the current filter view. </p> <p>  Deletes the current history filter view. </p>
Order	The sequence of actions that occurred.
Change Type	One of the predefined ChangeType values.
Context	Name of the specific value on which the action occurred.
Item Type	One of the predefined ItemType values.
Item Name	A user-defined name for the filter.
Old Value/New Value	Original parameter value and the changed value.
Details Key/Value	Identifying parameters and values for the selected history event.

Using Quick Acquisition

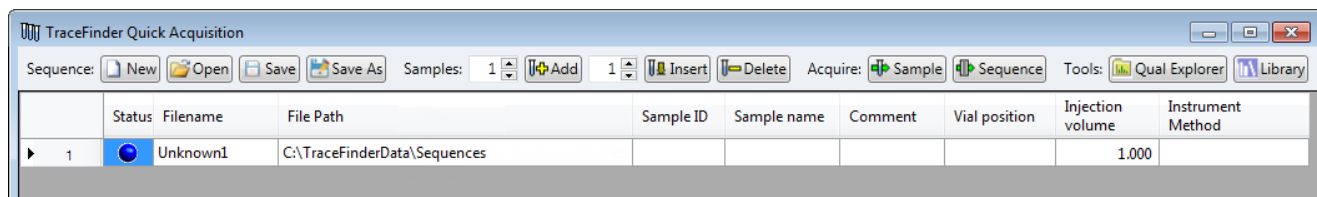
Use the quick acquisition feature to quickly submit samples from any mode in the application.

Note The Quick Acquisition feature is available only when you activate it in the Configuration console. Refer to Chapter 1, “Using the Configuration Console,” in the *TraceFinder Lab Director User Guide*.

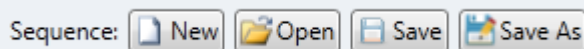
❖ To run a quick acquisition



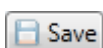
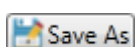
1. Choose **Tools > Quick Acquire Sample** from the main menu or click the **Quick Acquire Sample** icon, .

The TraceFinder Quick Acquisition window opens.

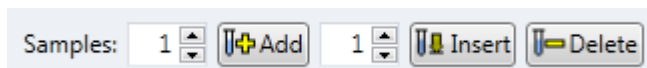





2. To create the sequence of samples that you want to acquire, do any of the following:
 - Use the Sequence buttons in the toolbar to open and save Xcalibur sequence (.sld) files.



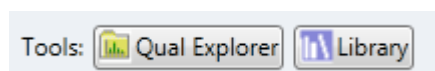
Button	Description
	Replaces the current sequence with a new sequence that contains one Unknown sample.
	Opens the Open dialog box where you can open a saved SLD file.
	Saves the current sequence as an SLD file in the C:\TraceFinderData\Sequences folder.
	Opens the Save As dialog box where you can save the current sequence to a new file name or location.

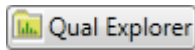

- Use the Samples buttons in the toolbar to create a sequence of samples.



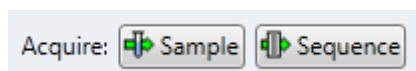
Button	Description
	Adds the specified number of new, empty samples to the end of the sample list.
	Inserts a new, empty sample or samples above the selected sample.
	Removes the selected samples from the sample list.

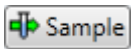
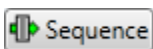
- (Optional) Use the Tools buttons in the toolbar to open a qualitative browser or the NIST library browser.



Button	Description
	Opens the qualitative explorer that you configured as your default qualitative explorer in the Configuration console. See Launching a Qualitative Explorer .
	Opens the NIST library browser. See Launching the NIST Library Browser .

- When you have completed your sequence of samples, click either of the Acquire buttons.



Button	Description
	Submits only the selected samples for acquisition, processing, or report generation.
	Submits the sequence for acquisition, processing, or report generation.

The application submits the samples for acquisition, processing, and report generation. See [Real Time Status – Acquisition Page](#).

❖ To see how to use the Quick Acquisition wizard

- Choose **Help > Animations**.
- From the list of animation topics, click **Quick Acquisition**.

Using Copy Down and Fill Down

This appendix describes the Copy Down and Fill Down commands that you can use to make entering column values easier.

- Use the Fill Down command for the Filename, Sample Name, Sample ID, and Vial Position columns.
- Use the Copy Down command for the Sample Type, Vial Position, Injection Volume, Conv Factor, Level, Comment, and other columns.

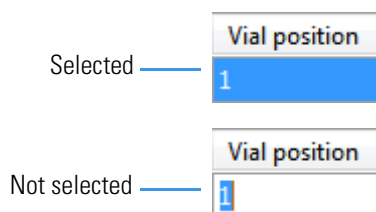
Follow these procedures:

- [To automatically copy column values](#)
- [To automatically enter sequential column values](#)
- [To use Copy Down or Fill Down for a range of samples](#)

❖ To automatically copy column values

1. Select the cell whose value you want to copy to all cells below it.

Observe the difference between a selected and nonselected cell.



2. Right-click and choose **Copy Down**.

The value is copied to all rows below the selected row.

❖ To automatically enter sequential column values

1. Enter a value for the first row of the fill down sequence.

This does not have to be the first sample row. You can begin the fill down procedure from any row in the sequence.

2. Select the cell whose value is the first in the fill down sequence.

Observe the difference between a selected and nonselected cell.



3. Right-click and choose **Fill Down**.

The application enters sequential column values starting with the value in the selected row and ending with the last row in the column.

You can repeatedly use the Fill Down command to create multiple sequences.

Vial position
A:A1
A:A2
A:A1
A:A2
A:A1
A:A2
A:A3
A:A4

When you use the Fill Down command for the Vial Position column with an autosampler configured, the application knows the number of vial positions configured in your autosampler and numbers the positions accordingly.

Vial position
A:A1
A:A2
A:A3
A:A4
A:A5
A:A6
A:B1
A:B2
A:B3
A:B4
A:B5
A:B6

❖ **To use Copy Down or Fill Down for a range of samples**

1. To select a range of sample values, do one of the following:

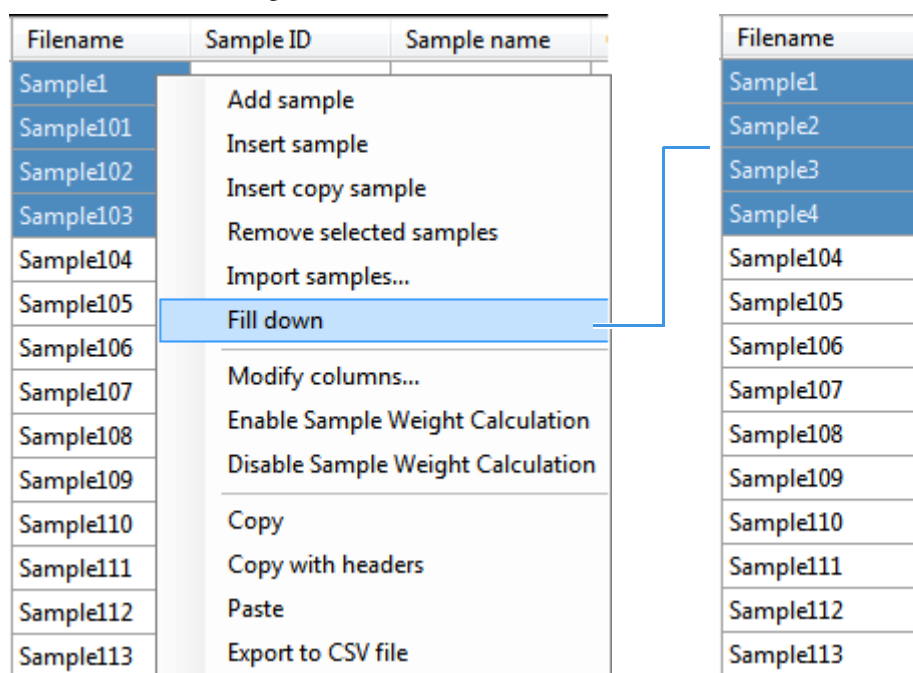
Drag your cursor to select a contiguous group of sample values.

—or—

Hold down the SHIFT key to select a contiguous group of sample values.

2. Right-click and choose the appropriate command from the shortcut menu.

The column values are copied or entered sequentially starting with the value in the first selected row and ending with the last selected row.



Moving Data Review Panes

This appendix describes the procedures you can use to move, dock, or float panes in all Data Review views. The procedures have accompanying animations.

Follow these procedures:

- [To move a docked pane](#)
- [To make a pane floating or dockable](#)
- [To change a pane from a docked pane to a tabbed pane](#)

❖ To move a docked pane

1. Grab the title bar of the pane and begin dragging the pane.

The application displays docking arrows.



2. Drag the pane over one of the arrows.

As you hold the cursor over a docking arrow, the application displays a blue region indicating where this arrow will place the pane.

3. Drop the pane onto one of the arrows.



This animation shows the various ways that you can use the docking mechanism to move a pane. To view the animation, click the filmstrip, and then right-click and choose **Full Screen Multimedia**. To stop the animation, press ESC.

❖ **To make a pane floating or dockable**

Do one of the following:

- To make a dockable pane floating, right-click the title bar of the pane and choose **Floating**.

While a pane is set as floating, you cannot use the docking arrows to dock it or make it a tabbed pane.

- To make a floating pane dockable, right-click the title bar of the pane and choose **Dockable**.



This animation shows how to switch a pane from docked to floating and back to docked. To view the animation, click the filmstrip, and then right-click and choose **Full Screen Multimedia**. To stop the animation, press ESC.

❖ **To change a pane from a docked pane to a tabbed pane**

1. Grab the title bar of the pane and begin dragging the pane.

The application displays docking arrows.



Drop the pane on the center location to create a tabbed pane.

2. Hold the cursor over the center of the docking arrows to display a blue region indicating the location of the tabbed pane.
3. Drop the pane over the center of the docking arrows.

Note To change a floating pane to a tabbed pane, you must first make the pane a dockable pane, and then you can make it a tabbed pane.



This animation shows how to change a pane from a docked pane to a tabbed pane and back to a docked pane. To view the animation, click the filmstrip, and then right-click and choose **Full Screen Multimedia**. To stop the animation, press ESC.