

Thermo

TraceFinder

Version 1.1

User Guide

XCALI-97357 Revision A

June 2011





© 2011 Thermo Fisher Scientific Inc. All rights reserved.

LCquan and TSQ Vantage are trademarks, and Xcalibur and TSQ Quantum 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.

Windows, Excel, and Windows Vista are registered trademarks of Microsoft Corporation in the United States and other countries. Adobe, Acrobat, Flash, and Reader are registered trademarks of Adobe Systems Inc. in the United States and other countries.

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.

Thermo Fisher Scientific Inc. makes no representations that this document is complete, accurate or errorfree and assumes no responsibility and will not be liable for any errors, omissions, damage or loss that might result from any use of this document, even if the information in the document is followed properly.

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, June 2011

Minimum software requirements: Thermo Xcalibur 2.1.0 SP1; Microsoft Windows XP Professional SP3; (optional) Thermo LC Devices 2.5

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

Contents

| | Preface | vii |
|-----------|--|------|
| | Related Documentation | |
| | Special Notices | viii |
| | System Requirements | viii |
| | System Activation | |
| | Contacting Us | |
| Chapter 1 | Introduction | 1 |
| | About the TraceFinder Application | 1 |
| | TraceFinder Summary of Features | 3 |
| | TraceFinder Workflow | 4 |
| | Reporting Features | 5 |
| | Standard Report Types | 5 |
| | Custom Report Types | 6 |
| | Target Screening Report Types | 6 |
| Chapter 2 | Getting Started | 7 |
| - | Installing the TraceFinder Application | 8 |
| | Installing the QED and NIST Libraries | 10 |
| | Launching the NIST Library Browser | 13 |
| | Launching the Qual Browser | 14 |
| | Converting Legacy Data | 15 |
| | Choosing a Mode | 22 |
| Chapter 3 | Using the Configuration Mode | 27 |
| | User Administration | 30 |
| | Choosing User Roles | 36 |
| | Project Administration | 39 |
| | Working with Drives | 40 |
| | Working with Projects | 42 |
| | Compound Datastore | 44 |
| | Opening and Saving a Datastore | 45 |
| | Adding Compounds, Quan Peaks, and Confirming Ions to a Datastore | 51 |
| | Choosing Experiment Types | 55 |

iv

| | Application Configuration | 56 |
|-----------|---|-----|
| | Specifying the Reports Configuration | 58 |
| | Specifying User Security | |
| | Enabling a Compound Datastore | 67 |
| | Specifying Configuration Defaults | |
| | Specifying Multiplexing Parameters | |
| | Specifying Detection Parameters | |
| | Enabling Target Screening | |
| Chapter 4 | Using the Method Development Mode | 85 |
| | Working with Master Methods | 88 |
| | Creating a New Master Method | 89 |
| | Editing a Master Method | 104 |
| | Creating a Method Template | 188 |
| | Importing Published Master Methods | 198 |
| | Exporting SRM Data | 199 |
| | Working with Instrument Methods | 200 |
| | Working with Development Batches | 207 |
| | Creating a Development Batch | 207 |
| | Editing Samples in a Development Batch | 210 |
| | Acquiring Samples in a Development Batch | 212 |
| | Viewing Raw Data Files in the Qual Browser | 213 |
| Chapter 5 | Using the Acquisition Mode | 215 |
| | Working with Batches | 216 |
| | Opening and Navigating the Acquisition Mode | 216 |
| | Creating Batches | 218 |
| | Real-time Display | 246 |
| | Acquisition Page | 247 |
| | Instrument Page | 248 |
| | Devices Page | 249 |
| | Queues Page | 253 |
| | Real-time Trace Display | 257 |
| | Sample Types | 259 |
| Chapter 6 | Using the Data Review Mode | 261 |
| | Working in the Batch View | 262 |
| | Batch View Panes | |
| | Creating a New Batch | |
| | Editing a Batch | 271 |
| | Editing Report Output Formats | |
| | Setting Compound Active Status | |
| | Submitting a Batch | 278 |

| | Working in Data Review View | 284 |
|------------|---|-------|
| | Samples List | . 285 |
| | Quan Mode | . 295 |
| | Qual Mode | . 313 |
| | Working in Report View | . 325 |
| | Viewing Reports | 326 |
| | Generating Reports | . 330 |
| | Working with Reports | . 333 |
| | Working with Active View | . 336 |
| | Working in the Local Method View | . 351 |
| A | Damanda | 252 |
| Appendix A | • | |
| | Specifying Reports | |
| | Standard Reports | |
| | Custom Reports | |
| | Target Screening Reports | |
| | Report Flags | |
| | Sample Standard Reports | |
| | Batch Report | |
| | Batch Report Rev 1 | |
| | Blank Report | |
| | Calibration Report. | |
| | Calibration Density 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 | |
| | High Density Internal Standard Report Long | |
| | High Density Sample Report 1 | |
| | High Density Sample Report 1 Long | |
| | High Density Sample Report 2 | |
| | High Density Sample Report 2 Long | |
| | High Density Sample Report 3 | |
| | High Density Sample Report 3 Long | |
| | Internal Standard Summary Report | |
| | Ion Ratio Failure Report | |
| | LCSLCSD Report (Environmental and Food Safety only) | |
| | Manual Integration Report | |
| | Method Detection Limit Report (Environmental and Food Safety only). | |
| | Method Report | |
| | Method Validation Report (Environmental and Food Safety only) | 392 |

Contents

vi

| | Index | 409 |
|------------|---|-----|
| | Enter Filter Criteria | 408 |
| Appendix C | Using Filter Criteria | 407 |
| Appendix B | Using Copy Down and Fill Down | 403 |
| | TIC Summary Report | 402 |
| | TIC Report | 401 |
| | Surrogate Recovery Report | 400 |
| | Solvent Blank Report | 399 |
| | Quantitation Report - 2 | 397 |
| | Quantitation Report | 396 |
| | MSMSD Report (Environmental and Food Safety only) | 395 |

Preface

The TraceFinder application is the newest application in the series of Thermo Scientific LC/MS analytical software.

Contents

- Related Documentation
- Special Notices
- System Requirements
- System Activation
- Contacting Us

To suggest changes to documentation or to Help

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



Related Documentation

The TraceFinder documentation includes Help that you can access from the application and also includes the *TraceFinder User Guide* and four quick reference guides as PDF files that you can access from the TraceFinder Help menu and the Windows Start menu.

❖ To view the TraceFinder user guide or quick reference guides

Go to **Start > All Programs > Thermo TraceFinder > Manuals** and choose one of the following documents:

- TraceFinder User Guide
- TraceFinder Administrator Quick Reference Guide
- TraceFinder Acquisition Quick Reference Guide
- TraceFinder Data Review Quick Reference Guide
- TraceFinder Shortcut Menus Quick Reference Guide
- TraceFinder Custom Reports Tutorial

❖ To open TraceFinder Help

From the TraceFinder window, choose **Help > TraceFinder Help.**

To find a particular topic, use the Help Contents, Index, or Search panes.

Special Notices

This guide includes the following types of special notices:

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.

System Requirements

Your system must meet these minimum requirements.

| System | Requirements |
|---|---|
| Computer | 2.33 GHz processor dual core with 2 GB RAM CD/R-ROM drive Video card and monitor capable of 1280 × 1024 resolution (XGA) 75 GB available on the C: drive NTFS format |
| Instruments (supported or required) | AS3000 autosampler |
| Software | Microsoft[™] Windows[™] XP Professional SP3 Microsoft Office 2007 SP2 or Excel[™] 2007 SP2 Microsoft .NET Framework 3.5 SP 1 Thermo Foundation[™] 1.1 (available on the Xcalibur 2.1.0 CD) Thermo Xcalibur[™] 2.1.0 SP1 Adobe[™] Reader[™] 9.0 NIST[™] 2008 |

System Activation

When you first start the TraceFinder application, a dialog box displays the number of days remaining on your 30-day free trial. If your free trial has expired, the License Activation window opens.



Note You can open the License Activation window at any time during your trial period by choosing **Help > License Activation** from the TraceFinder menu. If you already have a permanent license, a message tells you that your product is fully licensed.

Two types of licenses are available:

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

The evaluation version is full-featured and automatically expires 30 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 TraceFinder 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 questions regarding activation, contact Thermo Fisher Scientific Technical Support in San Jose, CA:

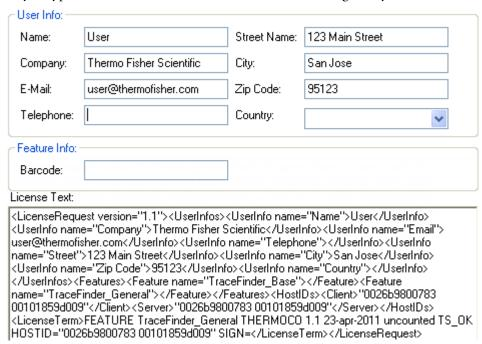
• E-mail: ThermoMSLicensing@thermofisher.com

• Fax: 408-965-6120

❖ To request an activation key

1. In the License Activation window, enter your information in the User Info area.

As you type, the License Text box creates an XML text string with your information.



2. In the barcode box, type the barcode printed on the TraceFinder CD.

The form of the barcode number is either xxxx-xxxx or xxxx-xxxx-xxxx-xxxx.

3. When you finish entering all your information, click **Copy**.

The application copies this XML text to the Clipboard.

If you have not completed all the information, a pop-up box opens, identifying the missing information.

4. Paste this XML text in the body of an e-mail and send the e-mail to ThermoMSLicensing@thermofisher.com.



To use your activation key

Note You must run the TraceFinder application with ITAdmin or LabDirector rights when entering the activation key.

- 1. When you receive your activation key, copy it from the e-mail.
- 2. Choose **Help > License Activation** from the TraceFinder menu.

The License Activation window opens.

3. Click Paste.

The application pastes the contents of the Clipboard to the License Text box.

4. Click Set.

The application is activated according to the type of authorization your license gives you.

Contacting Us

There are several ways to contact Thermo Fisher Scientific for the information you need.

❖ To contact Technical Support

Phone 800-532-4752 Fax 561-688-8736

E-mail us.techsupport.analyze@thermofisher.com

Knowledge base www.thermokb.com

Find software updates and utilities to download at mssupport.thermo.com.

To contact Customer Service for ordering information

Phone 800-532-4752 Fax 561-688-8731

E-mail us.customer-support.analyze@thermofisher.com

Web site www.thermo.com/ms

To get local contact information for sales or service

Go to www.thermoscientific.com/wps/portal/ts/contactus.

❖ To copy manuals from the Internet

Go to mssupport.thermo.com, agree to the Terms and Conditions, and then click **Customer Manuals** in the left margin of the window.

xii

❖ To suggest changes to documentation or to Help

- Fill out a reader survey online at www.surveymonkey.com/s/PQM6P62.
- Send an e-mail message to the Technical Publications Editor at techpubs-lcms@thermofisher.com.

Introduction

This chapter describes the general features of TraceFinder software.

Contents

- About the TraceFinder Application
- TraceFinder Summary of Features
- TraceFinder Workflow
- Reporting Features

About the TraceFinder Application

The TraceFinder application targets general quantitation and environmental markets, creating the workflows that laboratories use. It supports a focused workflow for specific non-bioanalytical laboratory use, instrument control, and method development functionality in a single software package. The TraceFinder application is the primary application for the TSQ Quantum™ XLS triple quadrupole mass spectrometers.

The TraceFinder application can export SRM data in .xml format so that other applications can import the files into their databases.

The TraceFinder application can import the following file types:

- Sample lists in .csv or .xml format
- Processing (.pmd) and instrument (.meth) method files from the Xcalibur data system
- Compounds from files that use the datastore (.xml) format
- Batches or methods from the TraceFinder 1.0.1 application

The TraceFinder 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 TraceFinder 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 software suite with specific instructions that enable scientific instruments to perform data acquisition.
- Processing method (.pmd): A proprietary file format for the Xcalibur software 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.

TraceFinder Directory Structure

The TraceFinder application creates folders for projects/subprojects/batches and templates in the ...\Thermo\TraceFinder directory. Within each batch folder, the application creates folders for data, methods, and reports.

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

Figure 1. Example batch directory structure



TraceFinder Summary of Features

The TraceFinder system provides a workflow-oriented approach to high-throughput quantitation. The system uses a batch-centric approach and tools to 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 both the general quantitation and the environmental and food safety applications. The application has a fully automated acquisition mode and a manual data review mode. You can use the data acquisition system to create and submit batches and monitor real-time review of results.

The TraceFinder 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 LabDirector, ITAdmin, Supervisor, Technician, and QAQC (quality assurance) roles
- Configuration mode for user administration, project administration, datastore administration, and application administration
- Method Development mode for editing instrument methods, setting processing and error flag parameters, and setting report options
- Acquisition mode that acts as a wizard to guide you in creating batches and running samples
- Data Review mode with batch views, data review, local method views, and report views
- Database-capable method development
- Quantification workflows, supporting capabilities present in the LCquan and LabForms applications
- Standard and customized report formats

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 TraceFinder application is structured with a typical laboratory workflow in mind. The user creates 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 lab. When you are ready to run a particular type of sample, select the appropriate method and begin.

When using the TraceFinder 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 how the raw data is acquired and processed, how the error checking information evaluates the results, and how the results appear in reports.

2. Create and submit a batch in the Acquisition mode.

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 the dashboard and 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 lower right corner of the TraceFinder window. This area displays a green, yellow, or red LED and a description of the number of samples in the queue (if any).

4. Evaluate the data in the Acquisition mode.

The Acquisition 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 Acquisition mode.

Use the Report View to view or print the reports for the currently selected 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 TraceFinder application can produce both standard reports and custom reports. The following types of reports meet the requirements of various methods and worldwide regulatory agencies, helping to track the performance of LC systems and methods. The reports divide into three groups: Standard, Custom, and Target Screening.

For additional information about standard, custom, and target screening reports and a sample of each standard report type, see "Reports" on page 353. Examples of standard reports (as PDFs) are also located in the C:\Thermo\TraceFinder\1.1\ExampleReports folder.

Standard Report Types

- Batch Report
- Batch Report Rev 1
- Blank 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
- High Density Internal Standard Report Long
- High Density Sample Report 1
- High Density Sample Report 1 Long
- High Density Sample Report 2
- High Density Sample Report 2 Long
- High Density Sample Report 3
- High Density Sample Report 3 Long
- Internal Standard Summary Report
- Ion Ratio Failure Report
- LCSLCSD Report (Environmental and Food Safety only)
- Manual Integration Report
- Method Detection Limit Report (Environmental and Food Safety only)
- Method Report
- Method Validation Report (Environmental and Food Safety only)
- MSMSD Report (Environmental and Food Safety only)
- Quantitation Report
- Quantitation Report 2
- Solvent Blank Report
- Surrogate Recovery Report
- TIC Report
- TIC Summary Report

1 Introduction Reporting Features

Custom Report Types

- AltCalibrationReport
- Alternate BatchReport
- Alternate CalibrationReport
- Alternate ConfirmationReport
- Alternate MatrixSpikeReport
- Alternate SampleReport
- Alternate SummaryReport
- BatchReport
- BlankReport
- CalibrationDensityReport
- CalibrationReport
- CheckStandardReport
- CompoundCalibrationReport
- ConfirmationReport
- ConfirmationReport2
- HighDensitySampleReport1Long
- HighDensitySampleReport2Long
- HighDensitySampleReport3Long
- HighDensitySampleReport4
- HighDensitySampleReport5
- QuantitationReport
- SteroidAnalysisReport

Target Screening Report Types

Target Screening reports are available only when you enable the Target Screening features. See "Enabling Target Screening" on page 83.

- Target Screening Long Report
- Target Screening Summary Report

7

Getting Started

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

Contents

- Installing the TraceFinder Application
- Installing the QED and NIST Libraries
- Launching the NIST Library Browser
- Launching the Qual Browser
- Converting Legacy Data
- Choosing a Mode

Installing the TraceFinder Application

Follow these instructions to install, start, and log on to the TraceFinder application.

To install the TraceFinder application

- 1. Follow the instructions on the included CDs to install the Xcalibur 2.1 data system and your instrument drivers.
- 2. Install the driver for your LC device and autosampler.
- 3. Insert the TraceFinder CD in the drive, and follow the on-screen instructions.

If the install windows do not automatically open, navigate to the Xinstall.exe file and launch it.

To start the TraceFinder application

1. Configure your instruments.



You cannot configure your instruments while the TraceFinder application is running.

2. Double-click the **TraceFinder** application icon on your desktop, or go to **Start > All Programs > Thermo TraceFinder > TraceFinder** or **TraceFinder EFS**.

By default, user security is not enabled and the application does not require a password. To enable user security, see "Specifying User Security" on page 66.

To log on to the TraceFinder application (when user security is enabled)

1. Enter your assigned user name.

Before you can log on to the TraceFinder application, a system administrator must set up a user account for you. The administrator assigns you a user name and password and gives you permission to access specific modes.

IMPORTANT If you are the administrator logging on for the first time with user security enabled, use **Administrator/Password** as the *username/password*.

2. Enter your password.

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



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

3. Click Login.

The TraceFinder dashboard opens. See "TraceFinder Dashboard" on page 25.

4. To exit the TraceFinder application without logging on, click Exit TraceFinder.

Figure 2. TraceFinder logon screen



Table 1. Logon screen parameters

| Parameter | Description |
|------------------|--|
| Username | The user's assigned user name. |
| Password | The assigned password for the user name. |
| Login | Verifies the user name and password, and displays the dashboard. |
| Exit TraceFinder | Closes the TraceFinder application without logging on. |

Installing the QED and NIST Libraries

When you are using triple quad instruments, follow these instructions to install the NIST and QED libraries.

❖ To install the NIST library

- 1. Launch the XInstall executable file.
- 2. Click the **NIST** button.

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 program folder, select **C:\Thermo**.

❖ To install the QED library

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

The Thermo Xcalibur Roadmap opens.



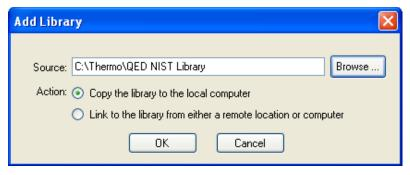
2. Select Tools > Library Manager.

The Thermo Library Manager dialog box opens. The NIST library is displayed in 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 application reports that it has added the library to the NIST software.

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

The application adds the QED library to the NIST Libraries list in the Library Manager dialog box.



2 Getting Started

Installing the QED and NIST Libraries

- 7. Click Exit in the Thermo Library Manager dialog box.
- 8. Start the TraceFinder application.
- 9. Go to the Method Development mode.
- 10. Click **Method View** in the navigation pane.
- 11. Choose **File > New > Method Template** from the main menu.

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



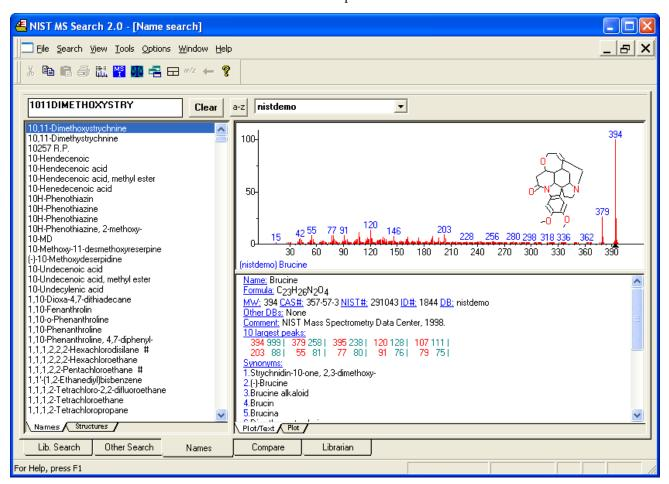
Launching the NIST Library Browser

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

❖ To open the NIST library browser

1. Choose **Go > Launch Library Browser** from the main menu.

The NIST MS Search window opens.



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

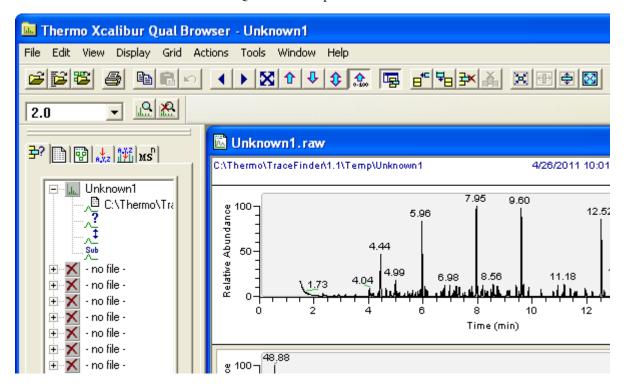
Launching the Qual Browser

Use the Qual Browser to view chromatograms and spectra from raw data files or qualitative processing results.

❖ To open the Qual Browser

1. Choose **Go > Launch Qual Browser** from the main menu.

The Thermo Xcalibur Qual Browser opens.



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

Converting Legacy Data

Use the TraceFinder Legacy Data Converter to convert methods, batches, method templates, or batch templates from a previous version of TraceFinder to the current version.

To open the TraceFinder Legacy Data Converter, choose **Go > Launch Legacy Data Converter** from the main menu.

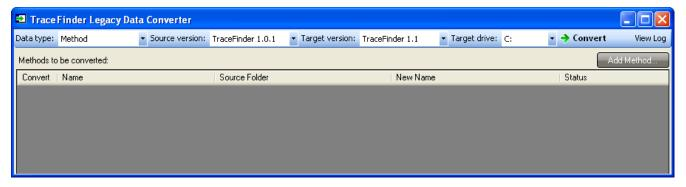
Use the following procedures:

- To convert a method
- To convert a batch
- To convert a method template
- To convert a batch template

❖ To convert a method

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

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



2. In the Source Version list, select a legacy version.

The legacy version defaults to TraceFinder 1.0.1.

- 3. In the Target Version list, select **TraceFinder 1.1**.
- 4. In the Target Drive list, select a fixed drive.

You cannot write the converted files to a mapped drive.

- 5. Click **Add Method** and browse to a method folder.
- 6. Select the method folder and click **OK**.

The application adds the method to the Methods to be Converted list:



2 Getting Started

Converting Legacy Data

- 7. Repeat step 6 for each method you want to convert.
- 8. (Optional) In the New Name column, type a new name for each method that you want converted.

If you leave this column blank, the converted method keeps the old name.

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

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 the method is successfully converted, the application writes the converted file to the C:\Thermo\TraceFinder\1.1\Methods folder.

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

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

===== Start Converting Method from TraceFinder 1.0.1 to TraceFinder 1.1 ========

Converting master method from 'C:\Thermo\TraceFinder\1.0.1\Method_4595\Method_4595.mmx'

Creating master method

Importing properties of object 'MethodData' from XML file 'C:\Thermo\TraceFinder\1.0.1\Method_4595\Meth Saving master method

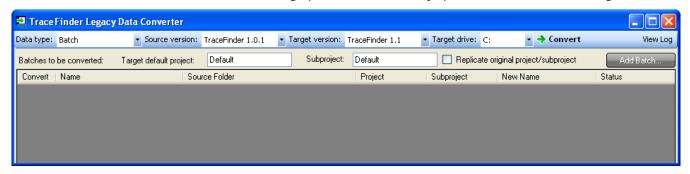
Copying instrument method

Successfully converted master method from 'C:\Thermo\TraceFinder\1.0.1\Method_4595\Method_4595.mmx' 'C:\Thermo\TraceFinder\1.1\Methods\Method_4595_new\Method_4595_new.mmx'

❖ To convert a batch

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

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



2. In the Source Version list, select a legacy version.

The legacy version defaults to TraceFinder 1.0.1.

3. In the Target Version list, select **TraceFinder 1.1**.

4. In the Target Drive list, select a fixed drive.

You cannot write the converted files to a mapped drive.

5. Do one of the following to create a project and subproject for the converted batch:

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.

- 6. Click Add Batch and browse to a batch folder.
- 7. Select the batch folder and click **OK**.

The application adds the batch to the Batches to be Converted list:



- 8. Repeat step 7 for each batch you want to convert.
- 9. (Optional) In the New Name column, type a new name for each batch that you want converted.

If you leave this column blank, the converted batch keeps the old name with "_1" appended.

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

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 the batch is successfully converted, the application writes the converted batch to the C:\Thermo\TraceFinder\1.1\Projects folder using either the original project and subproject names or the new names you entered.

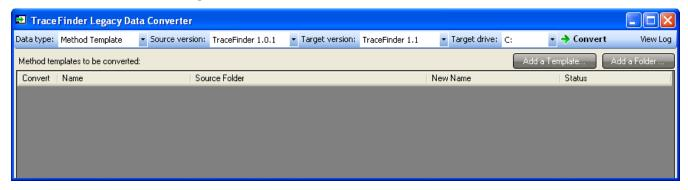
11. To view a log of the conversion, click **View Log**.

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

❖ To convert a method template

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

The TraceFinder Legacy Data Converter displays the interface for converting method templates.



2. In the Source Version list, select a legacy version.

The legacy version defaults to TraceFinder 1.0.1.

- 3. In the Target Version list, select **TraceFinder 1.1**.
- 4. In the Target Drive list, select a fixed drive.

You cannot write the converted files to a mapped drive.

- 5. To add a single method template, do the following:
 - a. Click **Add a Template** and browse to theC:\Thermo\TraceFinder\1.0.1\Templates\Methods folder.
 - b. Select the method template (.pmtx) file and click **Open**.

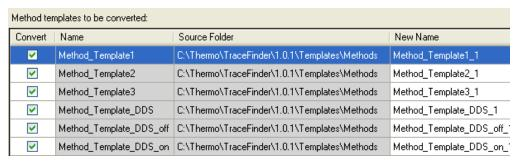
The application adds the method template to the Method Templates to be Converted list:



- 6. To add all the method templates in the Methods folder, do the following:
 - a. Click **Add a Folder** and browse to the C:\Thermo\TraceFinder\1.0.1\Templates folder.
 - b. Select the **Methods** folder and click **OK**.

Note The default folder for saved method templates is Templates\Methods. If you have templates in other folders, you can navigate to your folder.

The application adds all the method templates in the folder to the Method Templates to be Converted list:



- 7. Repeat step 6 for each method template you want to convert.
- 8. (Optional) In the New Name column, type a new name for each method template that you want converted.

If you leave this column blank, the converted template keeps the old name with "_1" appended.

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

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 C:\Thermo\TraceFinder\1.1\Templates\Methods folder.

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

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

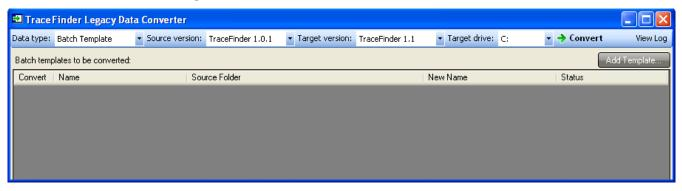
====== Start Converting Method Template from TraceFinder 1.0.1 to TraceFinder 1.1 ========= Converting method template from 'C:\Thermo\TraceFinder\1.0.1\Templates\Methods\Method_Template1.pmtx' Importing properties of object 'ProcMethodTemplateData' from XML file 'C:\Thermo\TraceFinder\1.0.1\Temp Saving the method template

Successfully converted method template from $C:\Thermo\TraceFinder\1.0.1\Templates\Method_Templates\Methods\Method_Template1_1.pmtx'$

To convert a batch template

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

The TraceFinder Legacy Data Converter displays the interface for converting batch templates.



2. In the Source Version list, select a legacy version.

The legacy version defaults to TraceFinder 1.0.1.

- 3. In the Target Version list, select **TraceFinder 1.1**.
- 4. In the Target Drive list, select a fixed drive.

You cannot write the converted files to a mapped drive.

- 5. Click **Add Template** and browse to the C:\Thermo\TraceFinder\1.0.1\Templates\Batches folder.
- 6. Select the batch template folder and click **OK**.

The application adds the batch template to the Batch Templates to be Converted list:



- 7. Repeat step 6 for each batch template you want to convert.
- 8. (Optional) In the New Name column, type a new name for each batch template that you want converted.

If you leave this column blank, the converted template keeps the old name with "_1" appended.

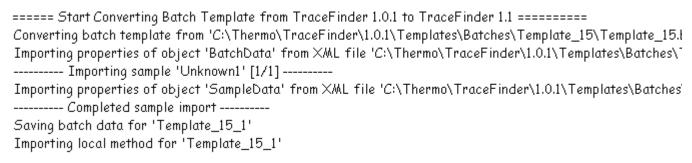
9. Select the check box for each batch template you will convert, and click **Convert**

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 C:\Thermo\TraceFinder\1.1\Templates\Batches folder.

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

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



Choosing a Mode

When user security is enabled, the dashboard provides the current user with options applicable to the role assigned to that user. The following table shows the available modes for each user role.

Table 2. User roles and permissions

| User role | Method Development | Acquisition | Data Review | Configuration |
|-------------|-----------------------|-------------|-------------|---------------|
| LabDirector | × | × | × | × |
| ITAdmin | | | | × |
| Supervisor | × | × | × | |
| Technician | | × | × | |
| QAQC | | | × | |

Note When user security is not enabled, all modes are available to all users.

Use the following procedures:

- To choose a mode
- To display a log of instrument errors
- To monitor the instrument status
- To watch the real-time display from the dashboard

❖ To choose a mode

1. From the dashboard, click the mode in which you want to work.

Your dashboard shows only the modes that you have permission to use. See "TraceFinder Dashboard" on page 25.

2. To change modes from within any of the TraceFinder application modes, click a mode button in the navigation pane.



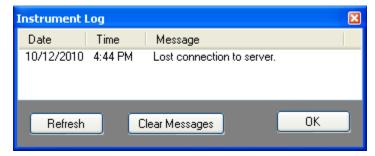
❖ To display a log of instrument errors

1. Right-click the status light in the lower right corner of any mode.



2. Select Instrument Log.

The Instrument Log dialog box opens.



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

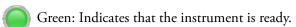
- 3. Do any of the following:
 - Click Refresh to display errors that occur after you open the Instrument Log dialog box.
 - Click Clear Messages to remove messages from the Instrument Log display.

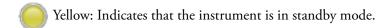
The application clears messages only from the Instrument Log display. These messages remain in the following log file:

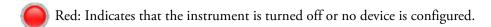
- C:\Thermo\TraceFinder\1.1\Logs\TraceFinder.log
- Click **OK** to dismiss the Instrument Log dialog box.

❖ To monitor the instrument status

To get immediate information about the instrument status, look at the LED in the lower right corner of the TraceFinder window.





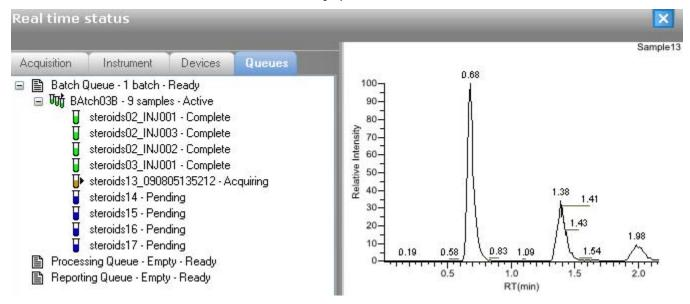


To watch the real-time display from the dashboard

Click Real Time Status.



The real-time status is displayed at the bottom of the dashboard.



For descriptions of all the features of the real-time display, see "Real-time Display" on page 246.

TraceFinder Dashboard

A dashboard for a user in the LabDirector role looks like this:

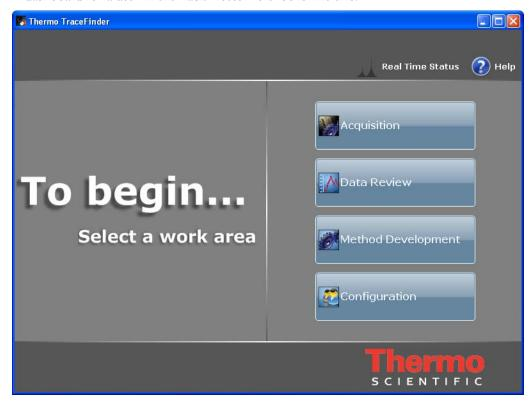


Table 3. TraceFinder dashboard screen parameters (Sheet 1 of 2)

| Parameter | Description |
|------------------|--|
| Real Time Status | Opens the real-time display for the current acquisition. The acquisition progress is displayed within the current mode window. |
| Help | Opens the TraceFinder Help. |
| Log Off | Logs off the current user and displays the logon screen. This function is available only when user security is enabled. See "Specifying User Security" on page 66. |
| Acquisition | Opens the Acquisition mode where you can create and review batches, batch data, reports, and local methods. See Chapter 5, "Using the Acquisition Mode." |
| Data Review | Opens the Data Review mode where you can review batches, batch data, reports, and local methods. See Chapter 6, "Using the Data Review Mode." |

2 Getting Started Choosing a Mode

Table 3. TraceFinder dashboard screen parameters (Sheet 2 of 2)

| Parameter | Description |
|--------------------|---|
| Method Development | Opens the Method Development mode where you can create a master method, an instrument method, or a development batch. See Chapter 4, "Using the Method Development Mode." |
| Configuration | Opens the Configuration mode where you can set permissions, assign users to roles, configure available reports and import new reports, and maintain the various databases, including the Compound Datastore. See Chapter 3, "Using the Configuration Mode." |

Using the Configuration Mode

This chapter discusses the configuration tasks assigned to the ITAdmin and LabDirector roles.

Contents

- User Administration
- Project Administration
- Compound Datastore
- Application Configuration

Users in the role of ITAdmin or LabDirector are responsible for the following:

- Handling the databases
- Applying roles to users
- Understanding security, users, and groups
- Creating local users and network groups
- Creating projects and subprojects
- Maintaining compounds in the compounds datastore

❖ To access the Configuration mode

Click **Configuration** from the dashboard or the navigation pane.



The Configuration mode navigation pane opens.

User Administration
Project Administration
Compound Datastore
Import Compounds
New Compound Datastore
Load Compound Datastore
Save Compound Datastore
Save As Compound Datastore
Recent files

Figure 3. Configuration mode navigation pane



Table 4. Configuration mode navigation pane functions (Sheet 1 of 2)

| Function | Description |
|------------------------|---|
| User Administration | Opens the User Administration view where you can add, remove, or edit user accounts and permissions. See "User Administration" on page 30. |
| | This task pane is available only when you have selected Enable User Security on the General page of the Application Configuration view. See "Application Configuration" on page 56. |
| Project Administration | Opens the Project Administration view where you can create and manage projects and subprojects. See "Project Administration" on page 39. |

Table 4. Configuration mode navigation pane functions (Sheet 2 of 2)

| Function | Description |
|---------------------------|--|
| Compound Datastore | Opens the Compound Datastore view where you can manage the definition of compounds in the current datastore. See "Compound Datastore" on page 44. |
| | This task pane is available only when you have selected Enable Compound Datastore on the General page of the Application Configuration view. See "Application Configuration" on page 56. |
| Application Configuration | Opens the Application Configuration view where you can specify available reports, user security, compound datastore, reporting defaults, multiplexing, detection algorithms, and target screening. See "Application Configuration" on page 56. |

User Administration

In the User Administration view of the Configuration mode, users in the role of LabDirector or ITAdmin can add, remove, or edit user accounts and permissions.

For detailed descriptions of each user role and the permissions and responsibilities for each role, see "Choosing User Roles" on page 36.

Use the following procedures:

- To open the User Administration view
- To add a user
- To edit user information
- To change a user password
- To remove a user

❖ To open the User Administration view

1. Click **Configuration** from the dashboard or the navigation pane.



The Configuration navigation pane opens.

Note The User Administration view is available only when you enable user security. Follow the instructions "To turn on user security" on page 66.

2. Click the **User Administration** task pane.



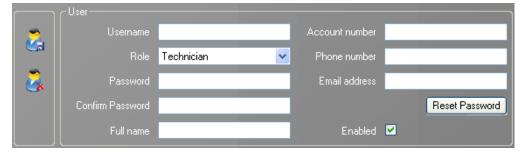
The User Administration view opens. See "User Administration view" on page 34.

To add a user

1. Click the **Add User** icon,



The application enables the parameters in the User area at the bottom of the view.



2. Type a unique name in the Username field.

3. Select a role from the Role list.

All users must be assigned to one of these defined roles. For detailed information about the permissions allowed for each role, see "Choosing User Roles" on page 36.

4. Type the user's password and type it again to the confirm it.

Make sure to communicate the password to the user.

- (Optional) Type the user's full name, account number, phone number, and e-mail address.
- 6. To enable this user logon, select the **Enabled** check box.

You can disable a user logon without deleting the user's information. Follow the instructions "To edit user information" on page 31.

7. Do one of the following:

When all the user information is correct, click the Save Changes icon,



The TraceFinder application adds the new user to the User Listing table, and the parameters in the User area are unavailable.

-Or-

To discard all information and not create a new user from the parameter values you entered, click the **Cancel Changes** icon,



The application discards all information and the parameters in the User area are unavailable.

❖ To edit user information

1. In the User Listing table, select a user.



Note Clicking anywhere in the row selects the user.

The user information populates the parameter fields in the User area.

User Administration

2. Click the Edit User icon,



The application enables the parameters in the User area.



3. Edit any of the parameter values.

If you are editing your own user name, the Enabled check box is unavailable because you cannot make your own account unavailable.

4. Do one of the following:



When all the user information is correct, click the Save Changes icon,

The TraceFinder application adds the new parameter values to the User Listing, and the parameters in the User area are unavailable.

- Or-



To discard all changes and not save the edits, click the Cancel Changes icon,

All changes are discarded, and the parameters in the User area are unavailable.

❖ To change a user password

1. In the User Listing table, select a user.

The user information populates the parameter fields in the User area.

2. Click the **Edit User** icon,



The parameters in the User area are enabled.



3. Click Reset Password.

The application makes the password and confirming password visible as a string of asterisks ******.

- 4. In the Password box, select the ***** and type a new password.
- 5. In the Confirm Password box, select the ***** and retype the new password.
- 6. Click the **Save Changes** icon, **?**

Make sure to communicate the new password to the user.

❖ To remove a user

1. In the User Listing table, select a user.

Note Clicking anywhere in the row selects the user.

The user information populates the parameter fields in the User area.

2. Click the **Remove User** icon,



If you select your current user name, the Remove User icon is unavailable. You cannot remove yourself.

3. When prompted, confirm that you want to remove this user.

If the user is currently logged on to the TraceFinder application, the user's current session is not affected.

4. Click OK.

Note Rather than completely removing the user, you can disable a user logon without removing all the user information from the system. Follow the instructions "To edit user information" on page 31.

3 Using the Configuration Mode

User Administration

Figure 4. User Administration view



Table 5. User Administration view parameters (Sheet 1 of 2)

| Parameter | Description | |
|-----------------|--|--|
| Security Groups | All permission levels defined in the TraceFinder application. For detailed descriptions of user permissions, see "Choosing User Roles" on page 36. | |
| User Listing | | |
| Username | User logon names. | |
| Role | The security group that defines user permissions. | |
| Account number | User account numbers. | |
| Phone number | User telephone numbers. | |
| Email address | User e-mail addresses. | |
| Enabled | Available or unavailable status for the user account. | |
| User | | |
| Username | Logon name for the current user. | |
| Role | Security group that defines the current user's permissions. | |
| Password | Logon password for the current user. | |
| Full name | The the current user's actual name. | |
| Account number | Optional account number for the current user. | |
| Phone number | Optional telephone number for the current user. | |

Table 5. User Administration view parameters (Sheet 2 of 2)

| Paramete | er | Description |
|-------------|----------------|--|
| Email ad | dress | E-mail address for the current user. Used to notify user of a randomly generated password. |
| Enabled | | Allows or disallows access for this user. When this user is currently logged on, disallowing takes effect after the user logs off. |
| Reset Pas | ssword | Makes the password visible as a string of asterisks so that you can select and change it. |
| lcon | Function | |
| | Add user | Enables the fields in the User area where you can enter information for a new user. |
| | Remove user | Deletes all information for the selected user. |
| > | Edit user | Enables the User area where you can edit any of the parameters for the selected user. |
| | Save changes | Adds the new parameter values to the User Listing table and disables the parameters in the User area. |
| | Cancel changes | Discards all new or edited information. |

Choosing User Roles

This section describes the responsibilities for five different user roles when user security is enabled: LabDirector, ITAdmin, Supervisor, Technician, and QAQC.

IMPORTANT User roles are in effect only when user security is enabled. When user security is not enabled, all users have access to all modes.

User Permissions

A laboratory director or an IT administrator assigns you to a role that gives you permission to access specific modes of the TraceFinder application. When you log on, the dashboard displays links to only the modes that you can access.

Table 6. User roles and permissions

| User role | Method Development | Acquisition | Data Review | Configuration |
|-------------|-----------------------|-------------|-------------|---------------|
| LabDirector | × | × | × | × |
| ITAdmin | | | | × |
| Supervisor | × | × | × | |
| Technician | | × | × | |
| QAQC | | | × | |

LabDirector

In the LabDirector role, you review graphically applicable data and manipulate data, batches, methods, and instruments.

A laboratory director is responsible for these tasks:

- Creating or editing methods for new levels of detection or adding new compounds to the existing database
- Reviewing data from the mass spectrometer
- Running samples and reviewing data collected by others
- Reporting the data
- Understanding the results and giving final approval of the released data before archiving

ITAdmin

In the ITAdmin role, you set security, manage users into roles, and manipulate the various databases. You are responsible for adding compounds into the various compound databases.

An IT administrator is responsible for these tasks:

- Handling the databases
- Applying roles to users
- Understanding security, users, and groups
- Creating local users and network groups

Supervisor

In the Supervisor role, you are responsible for setting up the instrument samples and using previously built sequences and methods for processing and acquiring data. You also develop and edit methods for processing and acquiring data, review the data, and distinguish between the need to rerun samples or pass reports up to the lab manager for final review. On a daily basis, you establish the priority for a list of samples to run and create the sequence of events.

A supervisor is responsible for these tasks:

- Submitting samples
- · Creating and submitting batches
- Reporting the data to management
- Creating or editing methods for new levels of detection or adding new compounds to the existing database
- Reviewing data from the mass spectrometer
- Understanding the results, who ran the batch, and who passed along the results before
 giving intermediate approval and sending the data to management
- Modifying new compounds or adjusting methods for specific result sets

Technician

In the Technician role, you are responsible for setting up the instrument samples and using previously built sequences and methods for processing and acquiring data. You also edit existing methods for processing and acquiring data, review collected data, and distinguish between the need to rerun samples or pass reports up to the supervisor. On a daily basis, you are responsible for gathering the list of samples to run and creating the sequence of events.

A technician is responsible for these tasks:

- Submitting samples
- Creating and submitting batches
- Creating data to be reviewed by management
- Receiving instructions for new sets of samples for the TraceFinder application to analyze after finishing the current analysis
- Reviewing data from the mass spectrometer
- Understanding the resulting data, making integration changes, and passing those changes up for further approval

3 Using the Configuration Mode

User Administration

QAQC

In the QAQC role, you review graphically applicable data and interpret the data, but you do not manipulate the data.

A QAQC technician is responsible for these tasks:

- Reviewing data from the mass spectrometer
- Understanding the results and who ran and passed along the results before giving intermediate approval and sending the data to management
- Receiving instructions for new sets of samples for the TraceFinder application to analyze after finishing the current analysis

Note In the QAQC role, you cannot submit batches for acquisition from the Acquisition mode.

Project Administration

In the LabDirector or ITAdmin role, you can create and manage projects and subprojects on fixed or network drives in the Project Administration view of the Configuration mode.

This section includes the following topics:

- Working with Drives
- Working with Projects
- ❖ To open the Project Administration view
 - 1. Click **Configuration** from the dashboard or the navigation pane.

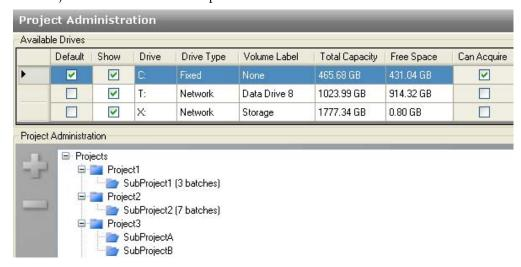


The Configuration navigation pane opens.

2. In the Configuration navigation pane, click Project Administration.



The Project Administration view opens.



By default, all projects are created under a main Projects folder on the C: drive:

C:\Thermo\TraceFinder\1.1\Projects

Working with Drives

Drives can be any of the following:

- Fixed: Directly connected to your computer.
- Network: Either a remote box or a mapped drive. A shared folder that is mapped to a drive letter might physically exist on your computer, but because it is mapped, it is considered to be a Network drive.
- Removable: Temporary drives such as a 3.5-inch disk or USB drive.

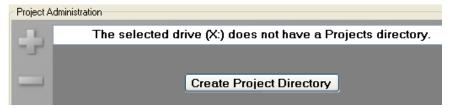
Use the following procedures:

- To choose a drive
- To change the default drive
- To hide a drive from display
- To refresh the drives display

❖ To choose a drive

1. In the Available Drives area, click any drive other than the default C: drive.

If you have not created a Projects directory on this drive, you see this message:



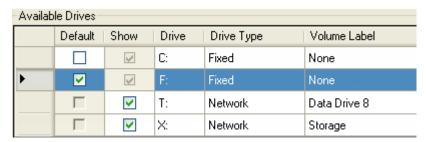
2. Click Create Projects Directory.

The TraceFinder application adds a new Projects directory to the selected drive. To create projects and subprojects on this drive, see "To create projects or subprojects" on page 42.

To change the default drive

Select the check box in the Default column.

Only fixed drives can be set as the default drive. The default drive is the first drive listed in all drive lists and is the only drive that you can use to acquire data.



❖ To hide a drive from display

Clear the check box in the Show column.

The application does not list the hidden drive in drive lists. You cannot hide the default drive.

To refresh the drives display

Right-click and choose Refresh from the shortcut menu.

The Available Drives table refreshes to show any drives that have changed, for example, if you have inserted a USB drive. You can now configure any new drives.

Working with Projects

When you create a batch, the application stores the data files, local method, and reports in a project and subproject that you create in the C:\Thermo\TraceFinder\1.1\Projects folder.

Use the following procedures:

- To create projects or subprojects
- To delete projects or subprojects
- To remove all empty folders
- To copy the folder hierarchy from another drive

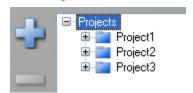
❖ To create projects or subprojects

1. Select the top-level project.

You can select the main Projects folder and create a new project under it, or you can select one of the existing projects and create a subproject under it.

When you select a project folder, the application enables the plus sign icon, indicating that you can create a folder within the selected folder.





2. Click the plus sign.

The TraceFinder application creates a new, unnamed project folder under the selected project.

3. While the new project is still highlighted, type a new name.

Project names are limited to 30 characters and can contain spaces and special characters, except for the following special characters: $\ \ / : *?" <> \ |$

Note After you add a subproject to a project, you cannot rename the project.

4. To save the new name, press ENTER or click anywhere in the view.

To delete projects or subprojects

1. Select the project or subproject you want to delete.



You can delete projects that do not have subprojects. You can delete subprojects that do not have batches. When the selected project or subproject is available for deletion, the application enables the minus sign icon,

- 2. Click the minus sign, or right-click and choose **Remove Project** from the shortcut menu.
- 3. At the prompt, click **Yes** to remove the selected project or subproject.

❖ To remove all empty folders

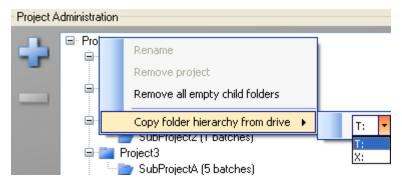
- 1. Select the project or subproject that contains empty folders.
- Right-click and choose Remove All Empty Child Folders from the shortcut menu.
 A dialog box asks if you want to remove all empty folders.

3. Click OK.

The application removes all folders that have no folders or files. There is no undo for this command.

❖ To copy the folder hierarchy from another drive

- 1. Select the top-level Projects directory in the Project Administration area.
- 2. Right-click and choose **Copy Folder Hierarchy from Drive** from the shortcut menu.
- 3. Choose a drive from the list of available drives.



At the prompt, you must confirm that you want to create a folder hierarchy that matches that of the specified drive.

4. Click OK.

To replicate the hierarchy from the specified drive, the application will add new folders to the current hierarchy, but it will not remove folders.

IMPORTANT The **Copy Folder Hierarchy from Drive** command copies only the project and subproject folders; it does not copy batches within the folders.

3 Using the Configuration Mode

Compound Datastore

Compound Datastore

In the LabDirector or ITAdmin role, you can manage compound definitions in the current datastore in the Compound Datastore view.

This section includes the following topics:

- Opening and Saving a Datastore
- Adding Compounds, Quan Peaks, and Confirming Ions to a Datastore
- Choosing Experiment Types

For a description of all the parameters in the Compound Datastore view, see "Compound Datastore View" on page 48.

Opening and Saving a Datastore

You can use the default datastore or you can create your own datastore.

Use the following procedures:

- To open the Compound Datastore editor
- To open a compound datastore
- To create a new compound datastore
- To save a datastore
- To save a datastore to a new name

❖ To open the Compound Datastore editor

1. Click **Configuration** from the dashboard or the navigation pane.

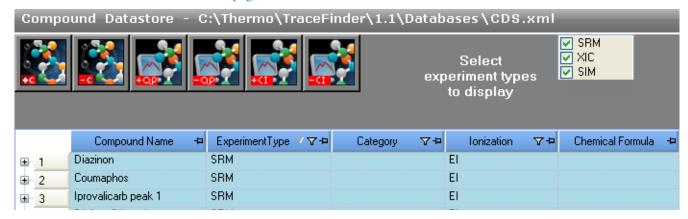


The Configuration navigation pane opens.

2. Click the **Compound Datastore** task pane.



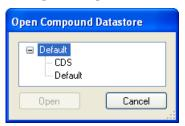
The current datastore opens in the Compound Datastore view. For a detailed list of all parameters and functions on the compound Datastore view, see "Compound Datastore View" on page 48.



To open a compound datastore

1. Click **Load Compound Datastore** in the Compound Datastore task pane.

The Open Compound Datastore dialog box opens.



2. Double-click the name of the datastore you want to open.

The selected datastore opens in the Compound Datastore view. See "Compound Datastore View" on page 48.

❖ To create a new compound datastore

Click **New Compound Datastore** in the Compound Datastore task pane.

A new, empty datastore opens in the Compound Datastore view.



You can import a file of compounds into the new datastore (following the instructions, To import compounds), or you can manually add compounds one at a time (following the instructions, To add a compound to the datastore).

❖ To save a datastore

1. Click **Save Compound Datastore** in the Compound Datastore task pane.

The application stores the database as

...\Thermo\TraceFinder\1.1\Databases\filename.xml

If the datastore contains any compounds that do not have associated quan peaks, the Invalid Compound Datastore Not Saved dialog box opens, listing the compounds.



2. To add a quan peak row to the listed compounds, click Continue.

The application returns you to the Compound Datastore view.

3. Add quan peaks to the incomplete compounds before you save the datastore.

See "To add a quan peak to a compound" on page 53.

❖ To save a datastore to a new name

1. Click **Save As Compound Datastore** in the Compound Datastore task pane.

The Save Compound Datastore dialog box opens.

- 2. Type a file name for the new compound datastore.
- 3. Click Save.

The application stores the database as

...\Thermo\TraceFinder\1.1\Databases\filename.xml

Figure 5. Save Compound Datastore dialog box

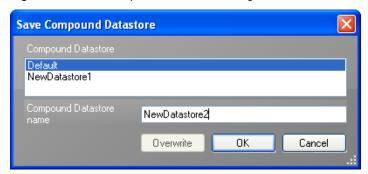


Table 7. Select Compound Datastore dialog box parameters

| Parameter | Description |
|--------------------|--|
| Compound Datastore | Lists the file name for the new datastore. |
| name | |
| Overwrite | Overwrites the selected datastore. |
| ОК | Writes the new datastore to the Databases folder. |
| Cancel | Closes the dialog box and makes no changes to the datastore. |

3 Using the Configuration Mode

Compound Datastore

Figure 6. Compound Datastore View

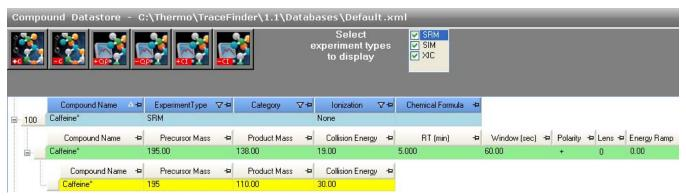


Table 8. Compound Datastore view parameters (Sheet 1 of 3)

| Parameter | Description | |
|--|--|--|
| Function icons | | |
| | Adds a new compound row. | |
| | Removes the selected row and all quan peak and confirming ion rows within it. | |
| ************************************** | Adds a new quan peak row to the compound. Each compound requires at least one quan peak. | |
| | Removes the selected row and all confirming ion rows within it. | |
| id. | Adds a new confirming ion row to the quan peak. | |
| -CT - | Removes the selected confirming ion row. | |
| Select experiment types | Specifies one of these experiment types that each use a different structure for the mass | |
| to display | filter. See "Choosing Experiment Types" on page 55. | |
| | SRM: Selected Reaction Monitoring | |
| | XIC: Extracted Ion Chromatogram | |
| | SIM: Single Ion Monitoring | |
| Compound parameters | | |
| Compound Name | Alphanumeric name assigned to the compound. | |
| Experiment Type | Experiment type: SRM, XIC, or SIM. | |
| Category | (Optional) Alphanumeric identifier. | |
| Ionization | (Optional) Alphanumeric identifier. | |
| | Valid values: None, ESI, APCI, EI, CI, APPI | |
| | Default: None | |
| Chemical Formula | (Optional) Alphanumeric chemical identifier. | |

Table 8. Compound Datastore view parameters (Sheet 2 of 3)

| Parameter | Description | |
|----------------------|--|--|
| Quan peak parameters | | |
| Precursor Mass | The mass-to-charge ratio of a precursor ion. The location of the center of a target precursor-ion peak in mass-to-charge ratio (<i>m</i> / <i>z</i>) units. In confirming ion rows, the precursor mass is a noneditable copy of the quan peak precursor mass. Available only for SRM experiments. Default: 0.0 Range: 10.000 to 2999.999 | |
| Product Mass | The mass-to-charge ratio of the quan ion. The location of the center of a target quan-ion peak in mass-to-charge ratio (m/z) units. Available only for SRM experiments. Default: 0.0 Range: 10.000 to 2999.999 | |
| Mass | The mass-to-charge ratio of the quan ion. The location of the center of a target quan-ion peak in mass-to-charge ratio (m/z) units. Available only for XIC and SIM experiments. Default: 0.0 Range: 10.000 to 2999.999 | |
| Collision Energy | The energy used when ions collide with the collision gas. Available only for SRM experiments. Range: -250.00 to 250.00 | |
| RT (min) | Retention time. The application uses RT and Window values to determine the start and stop time for the acquisition. Range: 0.00 to 999.00 Start time = RT – (Window/2) Stop time = RT + (Window/2) Start and stop range: 0.00 to 999.00 | |
| Window (sec) | The application uses RT and Window values to determine the start and stop time for the acquisition. Range: 0.00 to 499.50 Start time = RT – (Window/2) Stop time = RT + (Window/2) Start and stop range: 0.00 to 999.00 | |
| Polarity | + (positive) or – (negative) | |
| Lens | (Optional) Range: -400 to 400 | |
| Energy Ramp | (Optional) Available only for SRM experiments. Range: 0.00 to 200.00 | |

3 Using the Configuration Mode Compound Datastore

Table 8. Compound Datastore view parameters (Sheet 3 of 3)

| Parameter | Description | |
|--------------------------|---|--|
| Confirming ion parameter | s | |
| Precursor Mass | The mass-to-charge ratio of a precursor ion. The location of the center of a target precursor-ion peak in mass-to-charge ratio (m/z) units. Available as a noneditible field only for SRM experiments. Default: 0.0 Range: 10.000 to 2999.999 | |
| Product Mass | The mass-to-charge ratio of the quan ion. The location of the center of a target quan-ion peak in mass-to-charge ratio (m/z) units. Available only for SRM experiments. Default: 0.0 Range: 10.000 to 2999.999 | |
| Mass | The mass-to-charge ratio of the quan ion. The location of the center of a target quan-ion peak in mass-to-charge ratio (<i>m/z</i>) units. Available only for XIC and SIM experiments. Default: 0.0 Range: 10.000 to 2999.999 | |
| Collision Energy | The energy used when ions collide with the collision gas. Available only for SRM experiments. Range: -250.00 to 250.00 | |

Adding Compounds, Quan Peaks, and Confirming Ions to a Datastore

In the Compound Datastore view, you can import compounds into the datastore, add or remove compounds from the datastore, add quan peaks and confirming ions to a compound, or remove quan peaks or confirming ions from a compound.

Use the following procedures:

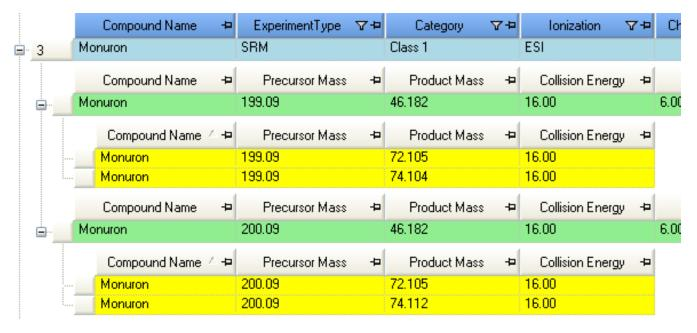
- To import compounds
- To add a compound to the datastore
- To remove a compound
- To add a quan peak to a compound
- To remove a quan peak
- To add a confirming ion to a quan peak
- To remove a confirming ion

❖ To import compounds

- 1. Click **Import Compounds** in the Compound Datastore task pane.
- 2. Browse to a .csv or .xml compounds file and click **Open**.

The TraceFinder application imports the compounds from the imported file, adds them to any compounds already in the datastore, and alphabetically sorts them.

When the application imports a compound that contains multiple quan peaks and confirming ions, it lists all the peaks under a single compound name, as in this example for Monuron:



❖ To add a compound to the datastore

1. Click the **Add Compound** button, or right-click the compounds list and choose **Add Compound** from the shortcut menu.

The application adds a new, empty compound row to the compounds table.



- 2. Click the first table cell, and enter the required Compound Name parameter.
- 3. (Optional) Change the value for the Experiment Type.

The default is SRM. For descriptions of available experiment types, see "Choosing Experiment Types" on page 55.

After you add a quan peak to the compound, you cannot change the experiment type, even if you cancel the quan peak.

4. (Optional) Type a value or select a value from the Category list.

You can use any alphanumeric string. After you type a new Category value, that value is available from the list.

5. (Optional) Change the values for Ionization.

The default is None.

6. (Optional) Type a value for the Chemical Formula.

You can use any alphanumeric string.

Each compound requires at least one quan peak.

To remove a compound

- 1. Select the compound row you want to delete.
- 2. Click the **Remove Compound** button, **Compound** from the shortcut menu.
- 3. If you are sure you want to delete the selected row, at the prompt, click **OK**.

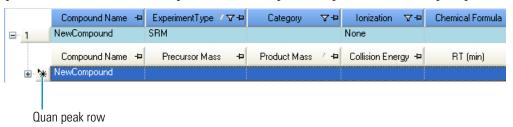
The application removes the selected row and all quan peak and confirming ion rows within it.

Tip If you add a row of compound information and do not complete all the column required values, you can right-click and choose **Cancel** to remove the entire row. You can cancel only an incomplete compound row.

To add a quan peak to a compound

- 1. Select the compound.
- 2. Click the **Add Quan Peak** button, with the shortcut menu.

The application adds a new quan peak row to the compound. A quan peak includes quantitative values for the compound. Each compound requires at least one quan peak.



3. Enter all required parameters.

For a list of required and optional parameters, see the list of "Compound Datastore view parameters" on page 48.

Tip You cannot add another new compound or save the compound datastore until you enter all required quan peak parameters or remove incomplete quan peaks from the compound.

4. Repeats steps 2 through 3 to add as many as six quan peaks to the compound.

❖ To remove a quan peak

- 1. Select the row of the quan peak you want to delete.
- 2. Click the **Remove Quan Peak** button, **Peak** from the shortcut menu.
- 3. If you are sure you want to delete the selected row, at the prompt, click **OK**.

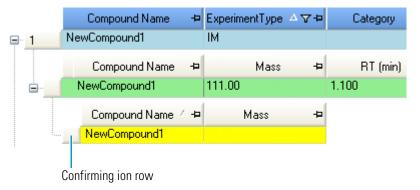
The application removes the selected row and all confirming ion rows within it.

Tip If you add a row of quan peak information and do not complete all the required values, you can right-click and choose **Cancel** to remove the entire row. You can cancel only incomplete quan peak rows.

To add a confirming ion to a quan peak

1. Click the **Add Confirming Ion** icon, or right-click the quan peak row and choose **Add Confirming Ion** from the shortcut menu.

The application adds a new confirming ion row to the quan peak. A confirming ion includes a mass value.



2. Type the required column values for the confirming ion.

The required confirming ion values differ for each experiment type. See "Choosing Experiment Types" on page 55.

3. Repeats steps 1 through 2 to add as many as 10 confirming ions to the quan peak.

To remove a confirming ion

- 1. Select the confirming ion row you want to delete.
- 2. Click the **Remove Confirming Ion** icon, confirming Ion from the shortcut menu.
- 3. If you are sure you want to delete the selected row, at the prompt, click **OK**.

The application removes the selected confirming ion row.

Tip If you add a row of confirming ion information and do not complete all the required values, you can right-click and choose **Cancel** to remove the entire row. You can cancel only incomplete confirming ion rows.

❖ To filter a list

Click the funnel icon, \mathbf{V} , in the column header.

For each column, the application displays filterable criteria (compound names, experiment types, categories, or ionization methods) in a list. In all lists, you can choose to view all criteria or a specific type of criterion for that column.

Choosing Experiment Types

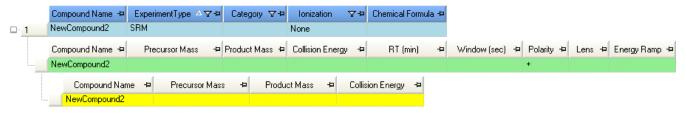
The TraceFinder application uses three experiment types: SRM, XIC, and SIM.

A compound datastore can include multiple experiment types for a single compound; however, each compound name and experiment type combination must be unique.

SRM – Selected Reaction Monitoring

The SRM experiment type supports triple quadrupole LC/MS. The mass filter includes precursor mass and narrow mass ranges to identify product masses. Imported compounds with no experiment type are treated as SRM data.

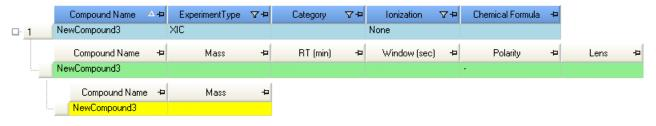
Confirming ions include values for product mass, collision energy, and a noneditable precursor mass.



XIC – Extracted Ion Chromatogram

The mass filter is a single, full scan which is post-processed to extract a peak for the ions of interest.

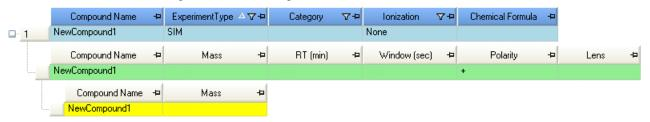
Confirming ions include a single mass value.



SIM - Single Ion Monitoring

The SIM experiment type supports single quadrupole LC/MS and Exactive[™] systems. The mass filter includes narrow mass ranges to identify product masses.

Confirming ions include a single mass value.



Application Configuration

In the LabDirector or ITAdmin role, you can enable features such as available reports, user security, compound datastore, reporting defaults, multiplexing, detection algorithms, and target screening. You can also choose the reports that are available to users, the application defaults, and the defaults used for peak detection.

This section includes the following tasks:

- Specifying the Reports Configuration
- Specifying User Security
- Enabling a Compound Datastore
- Specifying Configuration Defaults
- Specifying Multiplexing Parameters
- Specifying Detection Parameters
- Enabling Target Screening

To open the Application Configuration view

1. Click **Configuration** from the dashboard or the navigation pane.



The Configuration navigation pane opens.

2. Click **Application Configuration**.



The Application Configuration view opens.



Table 9. Application Configuration pages (Sheet 1 of 2)

| Page | |
|--------------------|--|
| Reports | See "Specifying the Reports Configuration" on page 58. |
| User Security | See "Specifying User Security" on page 66. |
| Compound Datastore | See "Enabling a Compound Datastore" on page 67. |

Table 9. Application Configuration pages (Sheet 2 of 2)

| Page | |
|-----------|--|
| Defaults | See "Specifying Configuration Defaults" on page 68. |
| Multiplex | See "Specifying Multiplexing Parameters" on page 70. |
| Detection | See "Specifying Detection Parameters" on page 71. |

3. In the Application Configuration list, click the type of information you want to configure.

Specifying the Reports Configuration

In the LabDirector or ITAdmin role, you can configure a list of reports that are available to users when they generate reports from the Method Development, Data Review, and Acquisition modes. From the Reports page, you can configure the standard, custom, or target screening reports.

Example PDFs of report formats are located in the following folder:

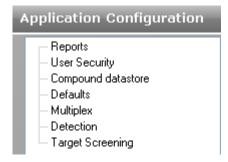
C:\Thermo\TraceFinder\1.1\ExampleReports

Use the following procedures:

- To open the Reports page
- To specify which reports are available
- To import new reports

To open the Reports page

In the Application Configuration view, click **Reports**.



The Reports page of the Application Configuration view opens.

- For a list of General Quantitation reports, see "General Quantitation Reports" on page 60.
- For a list of Environmental and Food Safety reports, see "Environmental and Food Safety Reports" on page 63.

To specify which reports are available

1. Use the directional arrows to move reports from the Installed Reports pane to the Displayed Reports pane.

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

In the Method Development, Data Review, and Acquisition modes, users can access all reports in the Displayed Reports pane.

2. To create a single composite report for an entire batch (rather than separate reports for each sample), select the **Batch Level** check box for the report.

Rather than creating separate reports for each sample, the application uses a composite of the data from all the samples to create a single report for the entire batch. Batch-level reports are prepended with a **B** to differentiate them.

Note Only reports with the ability to aggregate data at the batch level have the **Batch** Level check box enabled.

3. Do one of the following:

To apply the current selections, do the following:

a. Click **Apply**.

A message reminds you that you must restart the TraceFinder application before a user can access the reports you selected for the Method Development, Data Review, and Acquisition modes.

b. To restart the TraceFinder application now, click **Yes,** or to remain on the Reports page, click **No**.

-Or-

To return the report selections to their original state (when you first opened this page), click **Undo Changes**.

To import new reports

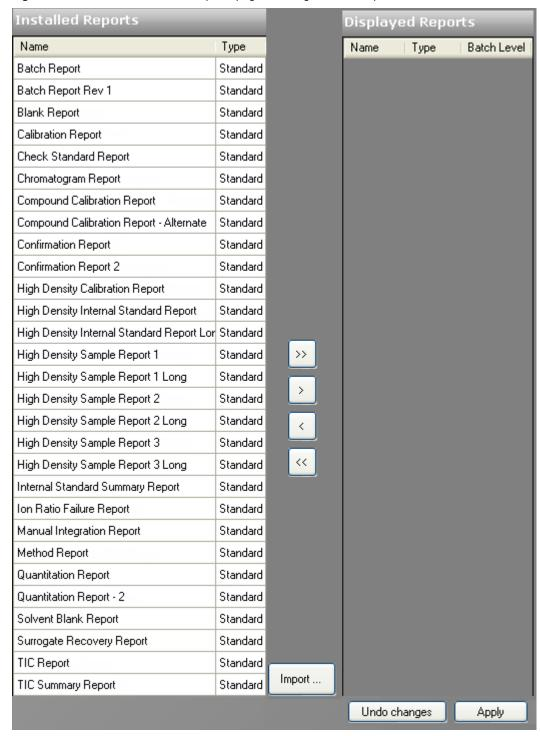
- 1. Click Import.
- 2. In the browser, locate a Crystal Reports .dll or Custom Reports .xltm file and open the file.

The application writes the imported report to the TraceFinder installation directory and displays the new report in the Installed Reports pane.

General Quantitation Reports

The General Quantitation configuration uses the following standard, custom, and target screening reports. For descriptions of the parameters on the Reports page, see "General Quantitation Reports page parameters" on page 62.

Figure 7. General Quantitation Reports page showing standard reports



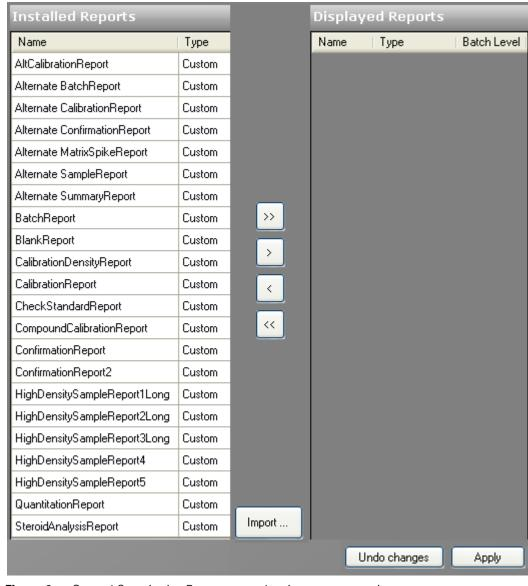
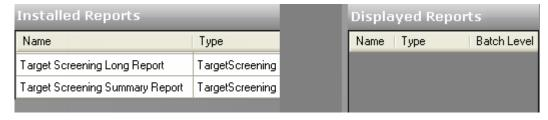


Figure 8. General Quantitation Reports page showing custom reports

Figure 9. General Quantitation Reports page showing target screening reports



Note Target Screening reports are available only when you enable target screening. See "Enabling Target Screening" on page 83.

3 Using the Configuration Mode

Application Configuration

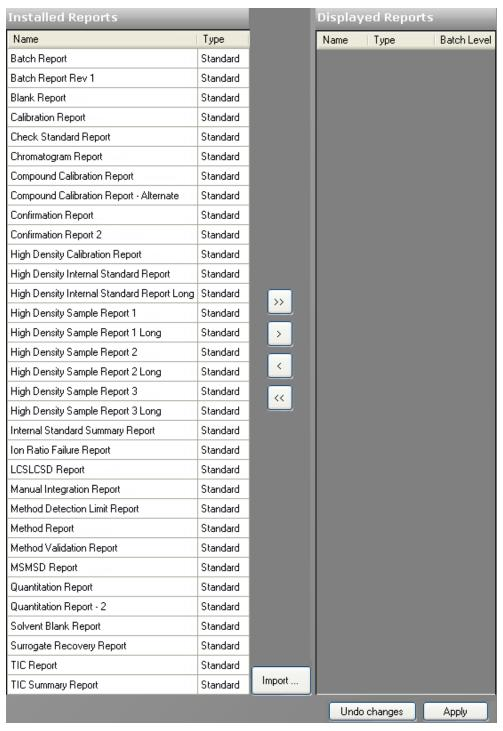
Table 10. General Quantitation Reports page parameters

| Parameter | Description |
|-------------------|---|
| Installed reports | All reports listed in the Installed Reports pane are potentially available but are not selected for use in the application. |
| Displayed reports | All reports listed in the Displayed Reports pane are selected for use in the application. |
| >> | Moves all reports from the Installed Reports list to the Displayed Reports list. |
| > | Moves the selected reports from the Installed Reports list to the Displayed Reports list. |
| < | Moves the selected reports from the Displayed Reports list to the Installed Reports list. |
| << | Moves all reports from the Displayed Reports list to the Installed Reports list. |
| Import | Opens a browser where you can select a report file to add to the Installed Reports list. |
| Undo changes | Returns the report selections to their original state (when you first opened this page). |
| Apply | Applies the current selections, and reminds you that you must restart the TraceFinder application before a user can access the reports you selected for the Method Development, Data Review, and Acquisition modes. |

Environmental and Food Safety Reports

The Environmental and Food Safety configuration uses the following standard, custom, and target screening reports. For descriptions of the parameters on the Reports page, see "Environmental and Food Safety Reports page parameters" on page 65.

Figure 10. Environmental and Food Safety Reports page showing standard reports



Installed Reports Displayed Reports Batch Level Name Type Name | Type AltCalibrationReport Custom Alternate BatchReport Custom Alternate CalibrationReport Custom Alternate ConfirmationReport Custom Alternate MatrixSpikeReport Custom Alternate SampleReport Custom Alternate SummaryReport Custom BatchReport Custom BlankReport Custom CalibrationDensityReport Custom CalibrationReport Custom CheckStandardReport Custom CompoundCalibrationReport Custom << ConfirmationReport Custom ConfirmationReport2 Custom HighDensitySampleReport1Long Custom HighDensitySampleReport2Long Custom HighDensitySampleReport3Long Custom HighDensitySampleReport4 Custom HighDensitySampleReport5 Custom QuantitationReport Custom Import ... SteroidAnalysisReport Custom Undo changes Apply

Figure 11. Environmental and Food Safety Reports page showing custom reports

Figure 12. Environmental and Food Safety Reports page showing target screening reports



Note Target Screening reports are available only when you enable target screening. See "Enabling Target Screening" on page 83.

Table 11. Environmental and Food Safety Reports page parameters

| Parameter | Description |
|-------------------|---|
| Installed reports | All reports listed in the Installed Reports pane are potentially available but are not selected for use in the application. |
| Displayed reports | All reports listed in the Displayed Reports pane are selected for use in the application. |
| >> | Moves all reports from the Installed Reports list to the Displayed Reports list. |
| > | Moves the selected reports from the Installed Reports list to the Displayed Reports list. |
| < | Moves the selected reports from the Displayed Reports list to the Installed Reports list. |
| << | Moves all reports from the Displayed Reports list to the Installed Reports list. |
| Import | Opens a browser where you can select a report file to add to the Installed Reports list. |
| Undo changes | Returns the report selections to their original state (when you first opened this page). |
| Apply | Applies the current selections, and reminds you that you must restart the TraceFinder application before a user can access the reports you selected for the Method Development, Data Review, and Acquisition modes. |

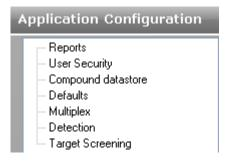
Specifying User Security

In the LabDirector or ITAdmin role, you can enable user security.

When user security is enabled, all users must log on to the TraceFinder application for access to only those modes assigned to their user role. See "Choosing User Roles" on page 36.

❖ To open the User Security page

In the Application Configuration view, click **User Security**.



The User Security page of the Application Configuration view opens.



To turn on user security

1. In the User Security area, select the **Enable User Security** check box.

When this check box is selected, all users must log on to the TraceFinder application for access to the modes assigned to their user role. See "Choosing User Roles" on page 36.

When this check box is cleared, users are not required to log on to the TraceFinder application. When they start the application, the dashboard is the first screen that users see and all modes are available to them. Only the User Administration view in Configuration mode is hidden and cannot be edited.

Note By default, user security is not enabled.

IMPORTANT If you are the administrator logging on with user security enabled, you can use **Administrator/Password** as the username/password.

2. In the Application Configuration view, click **Apply**.

A message reminds you that you must restart the TraceFinder application before the user security setting is enabled.

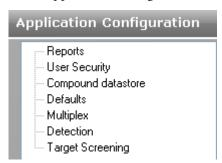
3. When you are prompted to restart the application, click **Yes**.

Enabling a Compound Datastore

In the LabDirector or ITAdmin role, you can enable the compound datastore.

To open the Compound Datastore page

1. In the Application Configuration view, click **Compound Datastore**.



The Compound Datastore page of the Application Configuration view opens.



2. In the Compound Datastore Selection area, select the **Enable Compound Datastore** check box.

Note By default, the Enable Compound Datastore option is not selected.

- 3. In the Application Configuration view, click **Apply**.
- 4. When prompted to restart the application, click **Yes**.

After restarting, the application implements the following changes:

- Displays the Acquisition List page on the Compounds page in the Master Method View. See "Editing the Compounds Page" on page 114.
- Displays the Compound Datastore task pane on the Configuration mode navigation pane. See "Compound Datastore" on page 44.
- Enables the Export SRM Data command in the Method Development mode. See "Exporting SRM Data" on page 199.

Specifying Configuration Defaults

In the Application Configuration view of the Configuration mode, as a user in the LabDirector or ITAdmin role, you can specify the default laboratory and instrument names, the mass precision, and the intensity scale to use for reporting.

Use the following procedures:

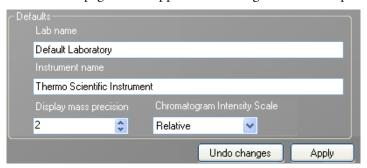
- To open the Defaults page
- To specify a default laboratory name and instrument name
- To specify default mass precision and the intensity scale

❖ To open the Defaults page

In the Application Configuration view, click **Defaults**.



The Defaults page of the Application Configuration view opens.



To specify a default laboratory name and instrument name

1. Type the name of your laboratory in the Lab Name box.

When you create a method, the application uses this default laboratory name for the Laboratory Name value on the General page of the Master Method View. The application uses this laboratory name in the report headings.

The application does not apply this default laboratory name to previously created methods. By default, the laboratory name is Default Laboratory.

2. Type the name of your instrument in the Instrument Name box.

When you create a batch, the application uses this default instrument name for the Instrument Name value. The application uses this instrument name in the report headings.

3. In the Application Configuration view, click **Apply**.

The application does not apply this default instrument name to previously created batches. By default, the instrument name is Thermo Scientific Instrument.

4. When prompted to restart the application, click **Yes**.

❖ To specify default mass precision and the intensity scale

1. In the Display Mass Precision box, set the mass precision decimal places value to an integer from 0 to 5, inclusive.

The default mass precision is 2. The TraceFinder application uses this mass precision value in the following locations:

- Reports:
 - Blank Report
 - Confirmation Report (data spectra, library spectra, quan ion display, and qual ion display)
 - All High Density reports (m/z values)
 - Ion Ratio Failure Report (quan ion and qual ion)
 - Manual Integration Report (m/z value)
 - Qualitative Summary Report (all m/z values)
 - Quantitation Report (QIon)
- All peaks on the Detection pages in the Method Development mode
- The spectrum display in the Data Review mode
- The spectrum display in the Method Forge dialog box
- 2. Select **Relative** or **Absolute** from the Chromatogram Intensity Scale list.

This sets the default display type for both quan and qual chromatograms displayed in data review and reports.

- 3. In the Application Configuration view, click **Apply**.
- 4. When prompted to restart the application, click **Yes**.

Specifying Multiplexing Parameters

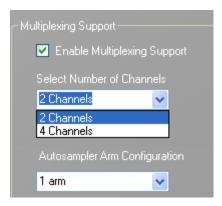
In the LabDirector or ITAdmin role, you can enable multiplexing and specify the maximum number of channels to use.

A user in the LabDirector, Supervisor, or Technician role can choose which channels to use for acquisition before submitting a batch from the Acquisition mode.

❖ To enable multiplexing

1. In the Application Configuration view, click **Multiplex**.

The Multiplexing Support page opens.



- 2. Select the **Enable Multiplexing Support** check box.
- 3. In the Select Number of Channels list, select 2 Channels or 4 Channels.
- 4. When you select 2 channels, select a 1- or 2-arm autosampler configuration.

The **1 arm** configuration enables channels 1 and 3; the **2 arm** configuration enables channels 1 and 2.

When you use a 4-arm configuration, autosampler 1 uses channels 1 and 3 and autosampler 2 uses channels 2 and 4.

- 5. To apply the multiplexing settings, do the following:
 - a. In the Application Configuration view, click **Apply**.
 - A message reminds you that you must restart the TraceFinder application before the multiplexing selections are reflected in the Acquisition mode.
 - b. To restart the TraceFinder application now, click **Yes,** or to remain on the Multiplexing Support page, click **No**.

Specifying Detection Parameters

In the LabDirector or ITAdmin role, you can specify detection parameters for Genesis, ICIS, or Avalon detection algorithms. Use the Peak Detection Defaults page to specify a peak detection algorithm and its options and to determine the area under a curve.

This section includes procedures for specifying the following detection algorithms:

- Genesis Detection Method
- ICIS Detection Method
- Avalon Detection Method

❖ To specify common detection parameters

1. In the Application Configuration view, click **Detection**.

The Peak Detection Defaults page opens. See "Common peak detection areas" on page 72.

- 2. In the Detector Type list, select a detector type.
- 3. In the Mass Tolerance area, do the following:
 - a. Select the units of measure you want to use.
 - b. In the Value box, specify the number of millimass units or parts per million to use as the upper limit.

The application applies this mass tolerance to the extracted chromatograms. The default is 500 MMU.

- 4. In the Retention Time area, do the following:
 - a. In the Window box, specify the width of the window (in seconds) to indicate how far around the expected retention time the system will look for a peak apex.
 - b. In the View Width box, specify the viewable size (in minutes) of the ion chromatogram display.
- 5. In the Ion Ratio Parameters area, do the following:
 - a. In the Window Type box, select **Absolute** or **Relative** as the calculation approach for determining the acceptable ion ratio range.
 - b. In the Window (+/-%) box, select the acceptable ion ratio range.
 - c. In the Ion Coelution box, select the maximum difference in retention time between a confirming ion peak and the quantification ion peak.
- 6. In the Peak Detection Parameters area, select one of the detection algorithms: **Genesis**, **ICIS**, or **Avalon**.

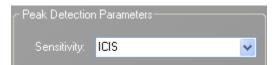


Figure 13. Common peak detection areas

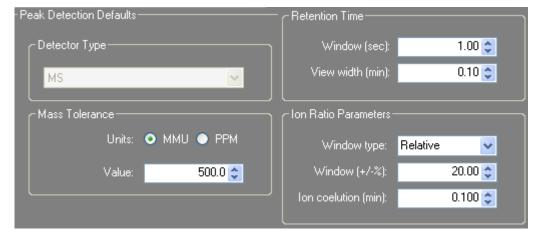


Table 12. Common peak detection parameters

| Parameter | Description |
|------------------------|---|
| Detector Type | |
| Mass Tolerance | |
| Units | (Default) MMU (millimass units) MMU is a static calculation to the extracted mass. PPM (parts per million) PPM is a variable calculation dependent on the actual mass. The smaller the mass, the narrower the tolerance range. The larger the mass, the wider the tolerance range. |
| Value | Upper limit of MMU or PPM. Default: 500 Range: 0.1 through 50 000 |
| Retention Time | |
| Window (sec) | Width of the window (in seconds) to indicate how far around the expected retention time the system will look for a peak apex. |
| View width (min) | Viewable size (in minutes) of the ion chromatogram display. Changing the view width does not affect the process of peak detection; the TraceFinder application uses it only for graphical display. |
| Ion Ratio Parameters | |
| Window type | The absolute or relative calculation approach for determining the acceptable ion ratio range. |
| Window (+/-%) | The acceptable ion ratio range. |
| Ion coelution (min) | The maximum difference in retention time between a confirming ion peak and the quantification ion peak. |

Genesis Detection Method

The TraceFinder application provides the Genesis peak detection algorithm for backward compatibility with Xcalibur 1.0 studies.

Figure 14. Genesis peak detection page

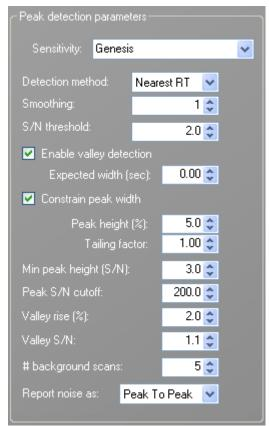


Table 13. Genesis peak detection page parameters (Sheet 1 of 3)

| Parameter | Description |
|------------------|---|
| Sensitivity | Specifies the Genesis peak detection algorithm. |
| Detection method | Highest peak: Uses the highest peak in the chromatogram for component identification. |
| | Nearest RT: Uses the peak with the nearest retention time in the chromatogram for component identification. |
| Smoothing | Determines the degree of data smoothing to be performed on the active component peak prior to peak detection and integration. The ICIS peak detection algorithm uses this value. |
| | Range: Any odd integer from 1 through 15 points Default: 1 |

Table 13. Genesis peak detection page parameters (Sheet 2 of 3)

| Parameter | Description |
|-------------------------|---|
| S/N threshold | Current signal-to-noise threshold for peak integration. Peaks with signal-to-noise values less than this value are not integrated. Peaks with signal-to-noise values greater than this value are integrated. |
| | Range: 0.0 to 999.0 |
| Enable valley detection | Uses the valley detection approximation method to detect unresolved peaks. This method drops a vertical line from the apex of the valley between unresolved peaks to the baseline. The intersection of the vertical line and the baseline defines the end of the first peak and the beginning of the second peak. |
| Expected width (sec) | The expected peak width parameter (in seconds). This parameter controls the minimum width that a peak is expected to have if valley detection is enabled. |
| | With valley detection enabled, any valley points nearer than the <i>expected width</i> /2 to the top of the peak are ignored. If a valley point is found outside the expected peak width, the TraceFinder application terminates the peak at that point. The application always terminates a peak when the signal reaches the baseline, independent of the value set for the expected peak width. |
| | Range: 0.0 to 999.0 |
| Constrain peak width | Constrains the peak width of a component during peak integration of a chromatogram. You can then set values that control when peak integration is turned on and off by specifying a peak height threshold and a tailing factor. Selecting the Constrain Peak Width check box enables the Peak Height (%) and Tailing Factor options. |
| Peak height (%) | A signal must be above the baseline percentage of the total peak height (100%) before integration is turned on or off. This text box is active only when you select the Constrain Peak Width check box. |
| | Range: 0.0 to 100.0% |
| Tailing factor | A factor that controls how the TraceFinder application integrates the tail of a peak. This factor is the maximum ratio of the trailing edge to the leading side of a constrained peak. This text box is active only when you select the Constrain the Peak Width check box. |
| | Range: 0.5 through 9.0 |

Table 13. Genesis peak detection page parameters (Sheet 3 of 3)

| Parameter | Description |
|-----------------------|---|
| Min peak height (S/N) | For the valley detection approximation method to use the Nearest RT Peak Identification criteria, this peak signal-to-noise value must be equaled or exceeded. For component identification purposes, the TraceFinder application ignores all chromatogram peaks that have signal-to-noise values that are less than the S/N Threshold value. |
| | Range: 0.0 (all peaks) through 999.0 |
| Peak S/N cutoff | The peak edge is set to values below this signal-to-noise ratio. |
| | This test assumes it has found an edge of a peak when the baseline adjusted height of the edge is less than the ratio of the baseline adjusted apex height and the peak S/N cutoff ratio. |
| | When the S/N at the apex is 500 and the peak S/N cutoff value is 200, the TraceFinder application defines the right and left edges of the peak when the S/N reaches a value less than 200. |
| | Range: 50.0 to 10000.0 |
| Valley rise (%) | The peak trace can rise above the baseline by this percentage after passing through a minimum (before or after the peak). This criteria is useful for integrating peaks with long tails. |
| | This method drops a vertical line from the apex of the valley between unresolved peaks to the baseline. The intersection of the vertical line and the baseline defines the end of the first peak and the beginning of the second peak. |
| | When the trace exceeds rise percentage, the TraceFinder application applies valley detection peak integration criteria. This test is applied to both the left and right edges of the peak. |
| | Range: 0.1 to 500.0 |
| Valley S/N | Specifies a value to evaluate the valley bottom. Using this parameter ensures that the surrounding measurements are higher. |
| | Range: 1.0 to 100.0 Default: 2.0 |
| # background scans | Number of background scans performed by the TraceFinder application. |
| Report noise as | Determines if the noise used in calculating S/N values is calculated using an RMS calculation or peak-to-peak resolution threshold. Options are RMS or Peak to Peak. |
| | |

ICIS Detection Method

The ICIS peak detection algorithm is designed for MS data and has superior peak detection efficiency at low MS signal levels.

Figure 15. ICIS peak detection page

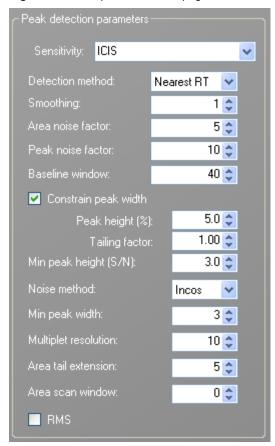


Table 14. ICIS peak detection page parameters (Sheet 1 of 3)

| Parameter | Description |
|------------------|---|
| Sensitivity | Specifies the ICIS peak detection algorithm. |
| Detection method | Highest peak: Uses the highest peak in the chromatogram for component identification. |
| | Nearest RT: Uses the peak with the nearest retention time in the chromatogram for component identification. |
| Smoothing | Determines the degree of data smoothing to be performed on the active component peak prior to peak detection and integration. The ICIS peak detection algorithm uses this value. |
| | Range: Any odd integer from 1 through 15 points Default: 1 |

Table 14. ICIS peak detection page parameters (Sheet 2 of 3)

| Parameter | Description |
|----------------------|--|
| Area noise factor | The noise level multiplier used to determine the peak edge after the location of the possible peak. The ICIS peak detection algorithm uses this value. |
| | Range: 1 through 500 Default: 5 |
| Peak noise factor | The noise level multiplier used to determine the potential peak signal threshold. The ICIS peak detection algorithm uses this value. |
| | Range: 1 through 1000 Default: 10 |
| Baseline window | The TraceFinder application looks for a local minima over this number of scans. The ICIS peak detection algorithm uses this value. |
| | Range: 1 through 500 Default: 40 |
| Constrain peak width | Constrains the peak width of a component during peak integration of a chromatogram. You can then set values that control when peak integration is turned on and off by specifying a peak height threshold and a tailing factor. Selecting the Constrain Peak Width check box enables the Peak Height (%) and Tailing Factor options. |
| Peak height (%) | A signal must be above the baseline percentage of the total peak height (100%) before integration is turned on or off. This text box is active only when you select the Constrain Peak Width check box. |
| | Range: 0.0 to 100.0% |
| Tailing factor | A factor that controls how the TraceFinder application integrates the tail of a peak. This factor is the maximum ratio of the trailing edge to the leading side of a constrained peak. This text box is active only when you select the Constrain the Peak Width check box. |
| | Range: 0.5 through 9.0 |

Table 14. ICIS peak detection page parameters (Sheet 3 of 3)

| Parameter | Description |
|-----------------------|---|
| Min peak height (S/N) | For the valley detection approximation method to use the Nearest RT Peak Identification criteria, this peak signal-to-noise value must be equaled or exceeded. For component identification purposes, the TraceFinder application ignores all chromatogram peaks that have signal-to-noise values that are less than the S/N Threshold value. |
| | Range: 0.0 (all peaks) through 999.0 |
| Noise method | The options are INCOS or Repetitive. |
| | INCOS: Uses a single pass algorithm to determine the noise level. The ICIS peak detection algorithm uses this value. |
| | Repetitive: Uses a multiple pass algorithm to determine the noise level. The ICIS peak detection algorithm uses this value. In general, this algorithm is more accurate in analyzing the noise than the INCOS Noise algorithm, but the analysis takes longer. |
| Min peak width | The minimum number of scans required in a peak. The ICIS peak detection algorithm uses this value. |
| | Range: 0 to 100 scans Default: 3 |
| Multiplet resolution | The minimum separation in scans between the apexes of two potential peaks. This is a criteria to determine if two peaks are resolved. The ICIS peak detection algorithm uses this value. |
| | Range: 1 to 500 scans Default: 10 |
| Area tail extension | The number of scans past the peak endpoint to use in averaging the intensity. The ICIS peak detection algorithm uses this value. |
| | Range: 0 to 100 scans Default: 5 |
| Area scan window | The number of allowable scans on each side of the peak apex. A zero value defines all scans (peak-start to peak-end) to be included in the area integration. |
| | Range: 0 to 100 scans Default: 0 |
| RMS | Specifies that the TraceFinder application calculate noise as RMS. By default, the application uses Peak To Peak for the noise calculation. RMS is automatically selected if you manually determine the noise region. |

Avalon Detection Method

The Avalon peak detection algorithm is designed for UV data. The Avalon peak detection algorithm also supports negative peaks.

Figure 16. Avalon peak detection page

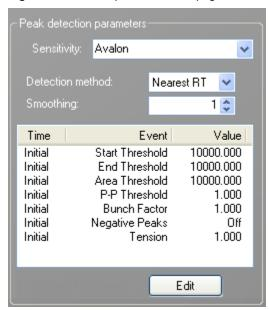


Table 15. Avalon peak detection page parameters

| Parameter | Description |
|------------------|---|
| Sensitivity | Specifies the Avalon peak detection algorithm. |
| Detection method | Highest peak: Uses the highest peak in the chromatogram for component identification. |
| | Nearest RT: Uses the peak with the nearest retention time in the chromatogram for component identification. |
| Smoothing | Determines the degree of data smoothing to be performed on the active component peak prior to peak detection and integration. The ICIS peak detection algorithm uses this value. |
| | Range: Any odd integer from 1 through 15 points Default: 1 |
| Time/Event/Value | Displays the events specified in the Avalon Event List dialog box. Initially displays only the default events that cannot be edited or deleted. |
| Edit | Opens the Avalon Event List dialog box where you can edit the Time/Event/Value parameters. See "Avalon Event List" on page 80. |

Avalon Event List

The event list includes both user-defined and noneditable default events. The application displays the default events when you choose Avalon sensitivity. You cannot delete these events or change their time or values. For a detailed list of events and value ranges, see "Event types" on page 81.

Figure 17. Avalon Event List dialog box

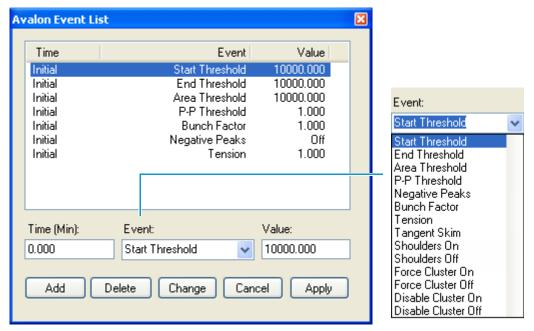


Table 16. Avalon Event List dialog box parameters

| Parameter | Description |
|------------|--|
| Time (Min) | Specifies the start time of the event. |
| Event | Specifies the type of event. For a detailed list of events and value ranges, see "Event types" on page 81. |
| Value | Specifies the value of the event. |
| Add | Adds a new event to the list with the current Time/Event/Value parameters. |
| Delete | Removes the selected Time/Event/Value parameter from the event list. |
| Change | Applies the current parameter values. |
| Cancel | Closes the dialog box without making any changes. Any additions, deletions, or changes revert to their original state. |
| Apply | Closes the dialog box. |

Figure 18. Event types



Table 17. Event type definitions (Sheet 1 of 2)

| Parameter | Description |
|-----------------|---|
| Start Threshold | Specifies the threshold at the start of a peak. The Start Threshold is directly related to the RMS noise in the chromatogram. Range: 0 to 999 999 999 |
| End Threshold | Specifies the threshold at the end of a peak. The End Threshold is directly related to the RMS noise in the chromatogram. Range: 0 to 999 999 999 |
| Area Threshold | Controls the area cutoff. Any peaks with a final area less than the area threshold will not be detected. This control is in units of area for the data. Range: 0 to 999 999 999 |
| P-P Threshold | The peak-to-peak resolution threshold controls how much peak overlap must be present before two or more adjacent peaks create a peak cluster. Peak clusters have a baseline drop instead of valley-to-valley baselines. Specified as a percent of peak height overlap. Range: 0.1 to 99.99 |
| Negative Peaks | Permits detection of a negative going peak. Automatically resets after finding a negative peak. Valid values: On or Off |
| Bunch Factor | Specifies the number of points grouped together during peak detection. This event controls the bunching of chromatographic points during integration and does not affect the final area calculation of the peak. A high bunch factor groups peaks into clusters. Range: 0 to 999 |

3 Using the Configuration Mode Application Configuration

Table 17. Event type definitions (Sheet 2 of 2)

| Parameter | Description | |
|---------------------|---|--|
| Tension | Controls how closely the baseline should follow the overall shape of the chromatogram. A lower tension traces the baseline to more closely follow changes in the chromatogram. A high baseline tension follows the baseline less closely, over longer time intervals. Range: 0 to 999.99 minutes | |
| Tangent Skim | Using this event, you can tangent skim any peak clusters. By default, it chooses the tallest peak in a cluster as the parent. You can also identify which peak in the cluster is the parent. Tangent skim peaks are detected on either side (or both sides) of the parent peak. Tangent skim automatically resets at the end of the peak cluster. Range: 0 to 1 | |
| Shoulders On | Allows peak shoulders to be detected (peaks which are separated by an inflection rather than a valley) Sets a threshold for the derivative. | |
| Shoulders Off | Disables peak shoulder detection. Range: 0 to 50 | |
| Force Cluster On | Force the following peaks to be treated as a cluster (single peak). | |
| Force Cluster Off | End the forced clustering of peaks. | |
| Disable Cluster On | Prevent any peaks from being clustered. | |
| Disable Cluster Off | Permit clusters to occur again. | |

Enabling Target Screening

In the Application Configuration view of the Configuration mode, you can enable target screening features.

Note You must have the ToxID application installed on your computer before you can generate Target Screening reports.

❖ To enable target screening

1. In the Application Configuration view, click **Target Screening**.

The Target Screening page of the Application Configuration view opens.



- 2. In the Target Screening area, select the **Enable Target Screening** check box.
- 3. In the Application Configuration view, click **Apply**.

A message reminds you that you must restart the TraceFinder application before the target screening feature is enabled.

4. When prompted to restart the application, click **Yes**.

Using the Method Development Mode

This chapter includes method development tasks assigned to the Supervisor or LabDirector roles.

Contents

- Working with Master Methods
- Working with Instrument Methods
- Working with Development Batches

From the Method Development mode, you can create a master method, an instrument method, or a simple development batch to test your instrument method.

❖ To access the Method Development mode

Click **Method Development** from the dashboard or the navigation pane.



The Method Development navigation pane opens. For a detailed description of all the features on the Method Development navigation pane, see "Method Development navigation pane" on page 86.



Figure 19. Method Development navigation pane

Table 18. Method Development navigation pane functions (Sheet 1 of 2)

| Function | Description | |
|---------------|---|--|
| Method View | See "Working with Master Methods" on page 88. | |
| Create method | Opens the Create Master Method dialog box where you can choose the process you want to use to begin your master method. | |
| Open method | Opens the Open Master Method dialog box where you can choose a master method to open. | |

Table 18. Method Development navigation pane functions (Sheet 2 of 2)

| Function | Description | | |
|----------------------------|---|--|--|
| Import Published Method | Opens the Import Published Method dialog box where you can select a published method to import. | | |
| Export SRM data | Writes the selected reaction monitoring (SRM) table to the following file: | | |
| | \Thermo\TraceFinder\1.1\Methods\methodname.xml | | |
| | You can use the data in this file in the instrument method editor when you open the TSQ 2.1 application. This command is available only when you have selected the Enable Compound Datastore option on the Compound Datastore page of the Application Configuration view. See "Application Configuration" on page 56. | | |
| | The compounds in the Acquisition List must contain at least one SRM experiment type. | | |
| Recent Files | Displays recently saved master methods. | | |
| Instrument View | See "Working with Instrument Methods" on page 200. | | |
| New instrument method | Opens the Instrument View where you can specify instrument settings for your configured instruments. | | |
| | If the instrument 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. | | |
| Open instrument method | Opens a browser where you can choose an instrument method to open. | | |
| Development Batch | See "Working with Development Batches" on page 207. | | |
| Select batch location | Specifies a location to store temporary development batch raw data files. | | |
| New sample list | Removes acquired samples from your development batch, so you can begin a new sample list. | | |
| Open Qual Browser | Opens the Thermo Xcalibur Qual Browser where you can view the acquired raw data files. | | |

Working with Master Methods

The TraceFinder application uses a master method to specify the nature and types of acquisition, processing, and reporting that occur with a batch of samples. When you are testing for compounds in an assay, you can create a method designed specifically for that type of application.

A master method contains a list of compounds and an initial set of information for detecting, processing, and reporting those compounds.

When you create a master method, the TraceFinder application uses the method to determine how the software works with a set of samples to provide a set of meaningful results. The application uses an instrument method to define how raw data is acquired. The rest of the master method defines how the raw data is processed, how the flags information evaluates the results, and how the reporting functionality defines the way your data and results are output into reports.

The TraceFinder application applies your master method to a batch, which is a list of one or more samples to be processed and reported. Together, the master method and batch provide a workflow-oriented approach to the data processing and information reporting for batches of samples.

To speed up the creation of master methods, you can create a method template. Using a method template helps you to develop methods faster because the TraceFinder application saves all of your commonly used method settings in a template, such as the number of confirming ions or the use of data-dependent data.

This section includes instructions for the following tasks:

- Creating a New Master Method
- Editing a Master Method
- Creating a Method Template
- Importing Published Master Methods
- Exporting SRM Data

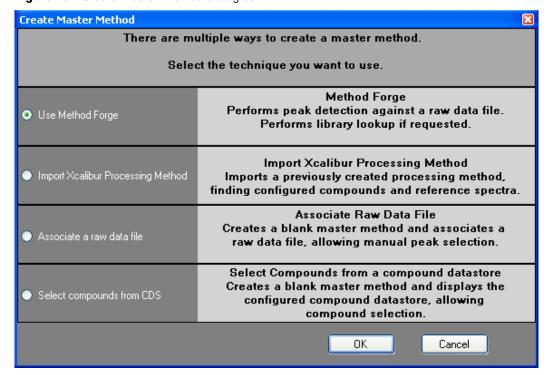
Creating a New Master Method

To begin a master method, follow any of four different procedures:

- Creating a New Method with Method Forge
- Importing an Xcalibur Master Method
- Associating a Raw Data File
- Selecting Compounds from the Compound Datastore

With each procedure, you begin the method in a specific way and then use the common features of the Master Method View to complete and save your master method.

Figure 20. Create Master Method dialog box



Creating a New Method with Method Forge

With Method Forge, you can create a new master method by manually selecting peaks, selecting multiple compounds, renaming peaks, or comparing mass spectra from the library searches. You can also choose to let the TraceFinder application automatically create a master method for you. For a detailed description of all the Method Forge parameters, see "Method Forge dialog box parameters" on page 96.

When the TraceFinder application automatically creates a master method for you, it performs the following functions:

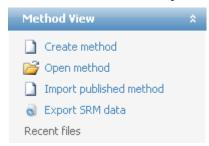
- Reviews your raw data file and identifies compounds that are present in your sample.
- Uses your mass spectral reference libraries to assign compound names and CAS numbers.
- Uses mass spectral information to select potential quantification and confirming ions and a reference mass spectrum for the compound.

Use the following procedures:

- To automatically select compounds to create a new method
- To manually select compounds to create a new method

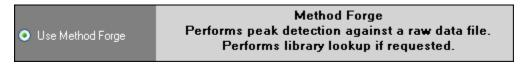
To automatically select compounds to create a new method

1. From the Method View task pane, click **Create Method**.



The Create Master Method dialog box opens.

2. Select the **Use Method Forge** option and click **OK**.



The Method Forge dialog box opens. For a detailed description of all the features on the Method Forge dialog box, see "Method Forge dialog box" on page 95.

Use the Method Forge to create a master method from an existing raw data file or to create a new raw data file to use for the master method.

3. In the Method Forge dialog box, do one of the following:

Select the **Use the Default Template** option.

-Or-

Select the **Select a Custom Template** option and highlight your custom template in the Method template table.

For detailed instructions on creating a custom method template, see "Creating a Method Template" on page 188.

- 4. Select the **Name the Master Method** check box and type a name for your master method. You can enter an existing method name and overwrite it when you create the method. If you do not specify a name, the application names the method for the raw data file used to create the method.
- 5. Select the **Automatically Create the Master Method** check box.
- 6. Do one of the following:
 - a. In the Raw File Selection area, choose **Use an Existing Raw Data File**.
 - b. Click the browse button and locate a raw data file to use for the method.
 - c. Go to step 8.

-Or-

- a. In the Raw File Selection area, choose **Acquire a New Raw Data File**.
- b. From the Instrument method list, select a method (.meth) file to use for acquiring the data.
- c. In the Raw Filename box, type the name of the file where the TraceFinder application will write the raw data file.
- d. In the Path box, type a path or click the browse button and locate a folder where the application will save the raw data file.
- e. (Optional) Type a comment about the acquired sample or the data file.
- 7. If you chose to acquire a new raw data file, do one of the following:

Choose Manual Injection.

-Or-

Do the following:

- a. Choose **Use Autosampler**.
- b. In the Vial Position box, type a vial position.
- c. In the Injection Volume box, type an injection volume.

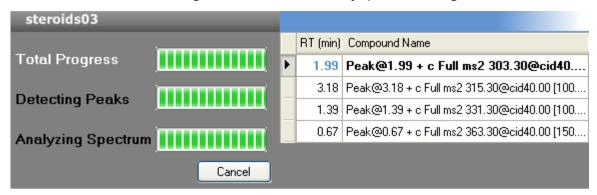
The minimum injection volume allowed is 0.1 $\mu L;$ the maximum injection volume allowed is 5000 $\mu L.$

4 Using the Method Development Mode

Working with Master Methods

8. To automatically create the master method, click \mathbf{OK} (or $\mathbf{Overwrite}$).

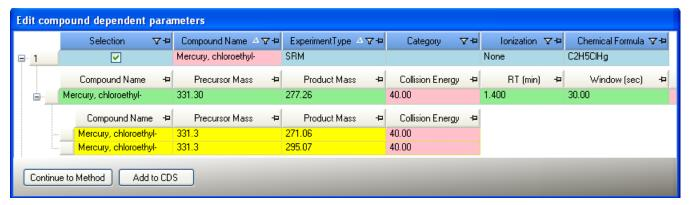
As the Method Forge creates the method, it displays the following status:



The Method Forge process searches the NIST library and displays the identified compound names instead of peak times.

When the acquisition completes, Method Forge performs peak detection, datastore searching, and characteristic ion and reference spectrum identification. Method Forge then loads this information into a new master method. This process occurs immediately if you selected a previously acquired raw data file.

If the compounds in the raw data file you used to create the method are not in the current compound datastore, the TraceFinder application displays the compounds in the Edit Compound Dependent Parameters dialog box.



9. (Optional) Select the compounds that you want to add to the compound datastore and click **Add to CDS**.

The selected compounds are added to the current compound datastore.

Note You must use the Add to CDS command before you continue to the method.

10. To use these compounds in your method and close the dialog box, click **Continue to Method**.

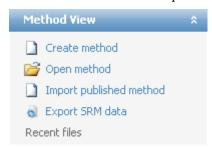
The TraceFinder application uses all compounds found in the raw data file in your method and displays the General page of the Master Method View. For a detailed description of all the features on the General page, see "General page" on page 111.

- 11. From the Instrument Method list on the General page, select an instrument method.
- 12. From the Qualitative Peak Processing Template list, select a method template for performing peak detection on quantitative samples following target compound analysis.
- 13. From the Background Subtraction Range Option list, select how you want the background subtraction range determined from one of these options:
 - **Before Peak**: Averages and subtracts a specified number of scans before the apex of the peak.
 - After Peak: Subtracts a specified number of scans following the apex of the peak.
 - **Both Sides of Peak**: Subtracts a specified number of scans from each side of the apex of the peak.
- 14. To save the new method, choose **File > Save** from the main menu.

For a detailed description of how to modify a master method, see "Editing a Master Method" on page 104.

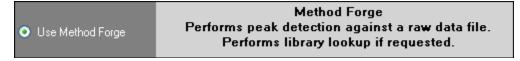
To manually select compounds to create a new method

1. From the Method View task pane, click Create Method.



The Create Master Method dialog box opens.

2. In the Create Master Method dialog box, select Use Method Forge and click OK.



The Method Forge dialog box opens. For a detailed description of all the features on the Method Forge dialog box, see "Importing an Xcalibur Master Method" on page 97.

3. In the Method Forge dialog box, do one of the following:

Select the **Use the Default Template** option.

−Or−

Select **Select a Custom Template** and highlight your custom template in the Method Template table.

For detailed instructions about creating a custom method template, see "Creating a Method Template" on page 188.

4 Using the Method Development Mode

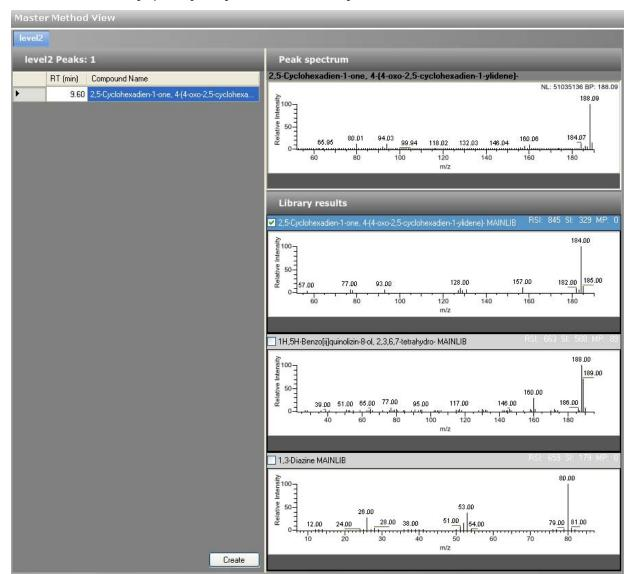
Working with Master Methods

4. Select the **Name the Master Method** check box and type a name for your master method.

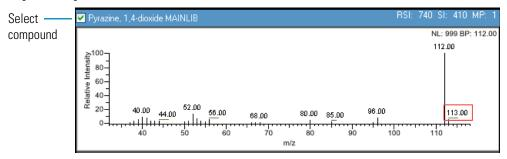
You can enter an existing method name and overwrite it when you create the method. If you do not specify a name, the method is named for the raw data file used to create the method.

- 5. Ensure that the **Automatically Create the Master Method** check box is not selected.
- 6. To select a raw data file, click the browse button and locate the file.
- 7. To manually create the master method, click **OK** (or **Overwrite**).

The Master Method View displays a list of possible matches in the Library Results pane. The TraceFinder application displays the best match in the Compound Name list and displays the peak spectrum for that compound.



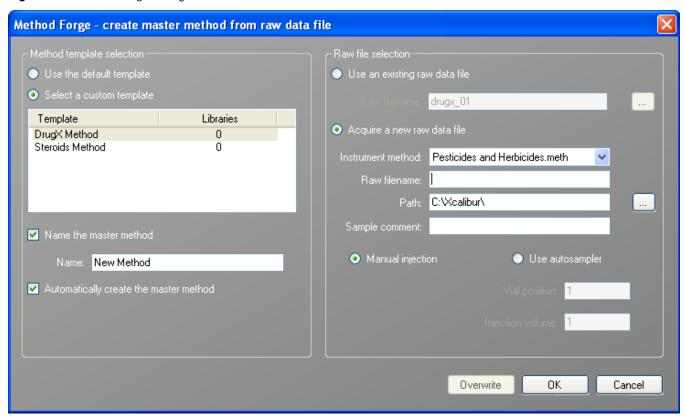
8. To use a compound other than the compound chosen by the TraceFinder application, scroll to the spectrum for that compound and select the compound name in the header of the spectrum pane.



- 9. After you manually select your compound, click **Create** to create the master method.
 - All compounds found in the raw data file are used in your method. The TraceFinder application displays the General page of the Master Method View. For a detailed description of all the features on the General page, see "General page" on page 111.
- 10. From the Instrument Method list on the General page, select an instrument method.
- 11. To save the new method, choose **File > Save** from the main menu.

For a detailed description of how to modify a master method, see "Editing a Master Method" on page 104.

Figure 21. Method Forge dialog box



4 Using the Method Development Mode

Working with Master Methods

Table 19. Method Forge dialog box parameters

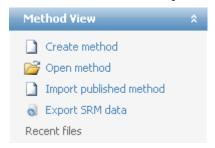
| Parameter | Description | | | |
|--|---|--|--|--|
| Method template selection | | | | |
| Use the default template | Creates a new method with the default template. | | | |
| Select a custom | Lists all the available method templates. | | | |
| template | For detailed instructions about creating a custom method template, see "Creating a Method Template" on page 188. | | | |
| Name the master method | The name for the new master method. | | | |
| Automatically create the master method | When the acquisition completes, Method Forge performs peak detection, library searching, and characteristic ion and reference spectrum identification. This information is loaded into a new master method. This process occurs immediately when you specify an existing raw data file. | | | |
| Raw file selection | | | | |
| Use an existing raw data file | Enables the Raw Filename box where you can select a raw data file to use to create the master method. | | | |
| Acquire a new raw data file | Enables functions to acquire data to create a raw data file to use to create the master method. | | | |
| Instrument method | The saved method (.meth) file to use for acquiring the data. | | | |
| Raw filename | The file name where the TraceFinder application writes the raw data. | | | |
| Path | The location where the TraceFinder application saves the raw data file. | | | |
| Sample comment | (Optional) Comment about the acquired sample or the data file. | | | |
| Manual injection | Performs a manual acquisition. | | | |
| Use autosampler | Performs an autosampled acquisition. | | | |
| Vial position | The tray vial number used for the autosampler acquisition. | | | |
| Injection amount | The volume (in milliliters) injected by the autosampler acquisition. | | | |
| Function button | | | | |
| Overwrite | Overwrites the specified master method name. This function is enabled only when the specified master method name already exists. | | | |
| OK | Creates a master method using the data and parameters you specified. | | | |
| Cancel | Closes the Method Forge and does not create a master method. | | | |

Importing an Xcalibur Master Method

You can create a new master method from an existing Xcalibur processing method.

❖ To import an Xcalibur master method

1. From the Method View task pane, click Create Method.

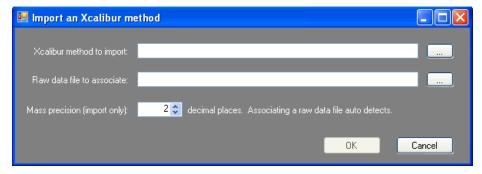


The Create Master Method dialog box opens.

2. In the Create Master Method dialog box, select the **Import Xcalibur Processing Method** option and click **OK**.



The Import an Xcalibur Method dialog box opens.



3. Click the browse button for the Xcalibur Method to Import box, browse to the Xcalibur processing method file, and open the file.

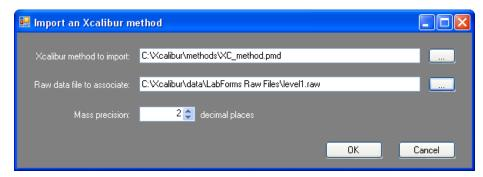
The TraceFinder application imports the compound information from the Xcalibur method file.

- 4. (Optional) Click the browse button for the Raw Data File to Associate box, browse to the raw data file, and open the file.
- 5. (Optional) Change the number of decimal places in the Mass Precision box.

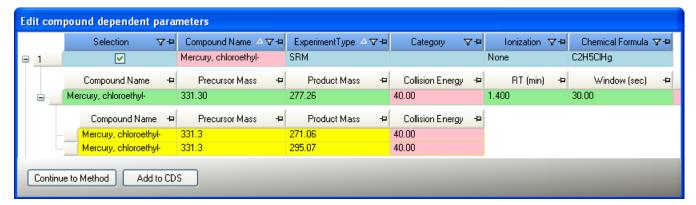
You can set the mass precision decimal places to any integer between 0 and 5, inclusive.

Note When you associate a raw data file, the application read the mass precision from the file and this parameter is unavailable.

6. Click OK.



If the compounds in the imported Xcalibur method file are not in the Compound Datastore, the TraceFinder application displays the compounds in the Edit Compound Dependent Parameters dialog box.



7. (Optional) Select the compounds you want to add to the Compound Datastore and click **Add to CDS**.

The selected compounds are added to the current compound datastore.

To add these compounds to the datastore, you must use the Add to CDS command before you continue to the method. When you click Continue to Method, the Edit Compound Dependent Parameters dialog box closes and you cannot return to add the compounds.

8. To add these compounds to your method and close the dialog box, click **Continue to Method**.

The TraceFinder application adds all compounds found in the imported Xcalibur method and displays the General page of the Master Method View. For a detailed description of all the features on the General page, see "General page" on page 111.

- 9. From the Instrument Method list on the General page, select an instrument method.
- 10. To save the new method, choose **File > Save** from the main menu.

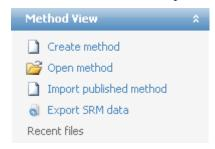
For a detailed description of how to modify a master method, see "Editing a Master Method" on page 104.

Associating a Raw Data File

You can use the compounds in a previously acquired raw data file to create a new master method.

Use the following procedures:

- To associate a raw data file with the method
- To add compounds to the method
- ❖ To associate a raw data file with the method
- 1. From the Method View task pane, click **Create Method**.

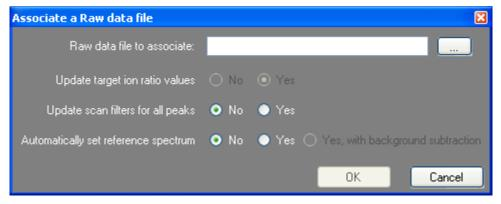


The Create Master Method dialog box opens.

2. In the Create Master Method dialog box, select the **Associate a Raw Data File** option and click **OK**.



The Associate a Raw Data File dialog box opens.



3. Click the browse button and locate a raw data file to associate with the method.

4 Using the Method Development Mode

Working with Master Methods

4. To set a reference spectrum, do one of the following:

Click Yes.

-Or-

Click Yes, with Background Subtraction.

This feature is available only when you have set background subtraction values on the General page of the Master Method View. See "Editing the General Page" on page 106.

The TraceFinder application adds all compounds found in the raw data file to your method and displays the General page of the Master Method View. For a detailed description of all the features on the General page, see "General page" on page 111.

- 5. From the Instrument Method list on the General page, select an instrument method.
- 6. To save the new method, choose **File > Save** from the main menu.

If the compounds in the associated raw data file are not found in the Compound Datastore, you cannot save the method. Follow the instructions "To add compounds to the method" on page 100.

For a detailed description of how to modify a master method, see "Editing a Master Method" on page 104.

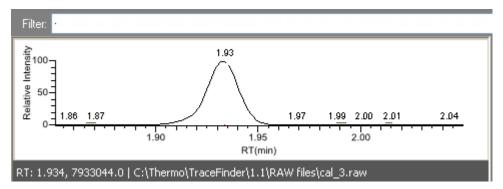
❖ To add compounds to the method

1. Click the **Compounds** tab.

The Detection page is selected by default.

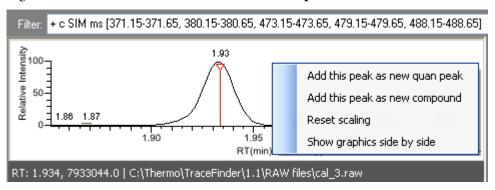


The Detection page shows an empty Compound list and displays the chromatographic data for the compounds in the raw data file.



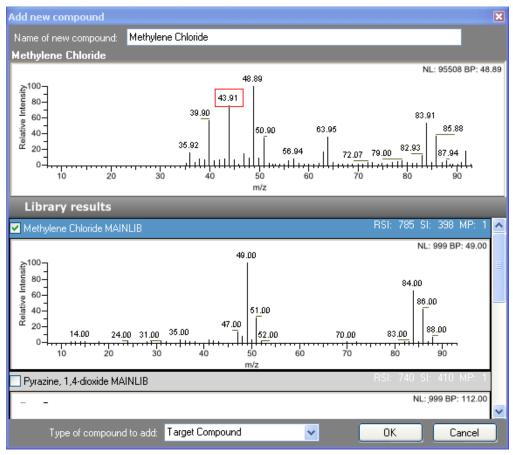
2. Select a filter from the Filter list.

- 3. Click the peak in the chromatogram that represents the compound you want to add to the method.
- 4. Right-click and choose **Add This Peak as New Compound** from the shortcut menu.



The TraceFinder application performs a library search for the selected compound. The application uses the first match it finds as the compound name, the base peak of the mass spectrum as the quan peak, and the second and third largest ions as the confirming ions.

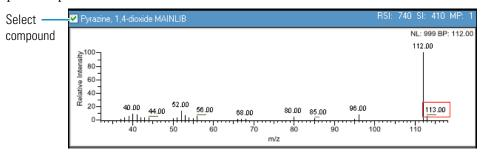
If the name of the first match is already in the library, the Add New Compound dialog box opens.



4 Using the Method Development Mode

Working with Master Methods

- 5. (Optional) In the Add New Compound dialog box, do the following:
 - a. To use a compound other than the compound already in the library, scroll to the spectrum for that compound and select the compound name in the title bar of the spectrum pane.



- b. In the Type of Compound to Add list, select a compound type.
- c. Click OK.
- 6. Repeat this procedure for each compound you want to add to the method.

For a detailed description of all the features on the Detection page, see "Editing the Compounds Page" on page 114.

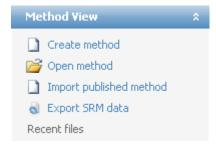
For a detailed description of how to modify a master method, see "Editing a Master Method" on page 104.

Selecting Compounds from the Compound Datastore

You can select compounds from the compound datastore to create a new master method. This method for creating a master method is available only when the compound datastore is enabled. See "Enabling a Compound Datastore" on page 67.

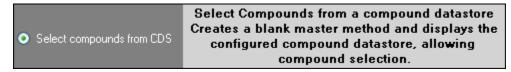
❖ To select compounds from the datastore

1. From the Method View task pane, click **Create Method**.

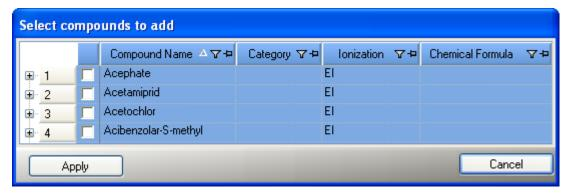


The Create Master Method dialog box opens.

In the Create Master Method dialog box, select the Select Compounds from CDS option and click OK.



The Select Compounds to Add dialog box opens, listing all the compounds defined in the compound datastore.



- 3. Select the check box for each of the compounds you want to add to the method.
- 4. To quickly select multiple compounds, right-click and use any of the commands on the shortcut menu.

Table 20. Select Compounds to Add shortcut menu

| Command | Description |
|--------------|--|
| Select All | Selects all compounds in the compound datastore. |
| Deselect All | Clears all compounds in the compound datastore. |
| Copy Down | Copies the state (checked or cleared) of the currently selected compound to all compounds below it in the compound list. |

5. Click Apply.

The TraceFinder application adds the selected compounds to the method.

Note After you add a compound to a method, the compound is no longer enabled in the Select Compounds to Add dialog box. You cannot remove the applied compounds from the method by returning to this dialog box. To remove a compound from a method, see "Acquisition List" on page 114.

- 6. From the Instrument Method list on the General page, select an instrument method.
- 7. To save the new method, choose **File > Save** from the main menu.

For a detailed description of how to modify a master method, see "Editing a Master Method" on page 104.

Editing a Master Method

You can open a master method to view or edit the compounds, method instructions, and reporting options in the method.

This section includes instructions for the following tasks:

- Opening a Master Method
- Editing the General Page
- Editing the Compounds Page
- Editing the Flags Page
- Editing the Groups Page
- Editing the Reports Page

Opening a Master Method

Use the TraceFinder application to open a master method that was created and saved in the current TraceFinder application or imported from TraceFinder version 1.0.1.

Use the following procedures:

- To open a saved master method
- To open a saved master method from the Recent Files list
- To open a saved master method
- 1. Click **Method Development** from the dashboard or the navigation pane.



The Method Development navigation pane opens.

In the Method View task pane, click Open Method, or choose
 File > Open > Master Method from the main menu.

The Open Master Method dialog box opens.



The Open Master Method dialog box displays all available methods. The list of TraceFinder methods can include any method created with TraceFinder 1.0.1 or greater.

3. Select a master method and click **Open**.

The General page for the selected method opens in the Method View. For a detailed description of all the features on the General page, see "General page" on page 111.

When you open a TraceFinder 1.0.1 method, the TraceFinder application copies all components of the selected method including its associated instrument method.

- If the TraceFinder 1.0.1 master method does not have an associated instrument method, the Master Method View for the imported method indicates that you must select an instrument method.
- If you open a TraceFinder 1.0.1 master method that uses an instrument method name already used in your Xcalibur\methods folder, you are prompted to overwrite the instrument method or use the method with the same name already in your methods folder.

To open a saved master method from the Recent Files list

1. Click **Method Development** from the dashboard or the navigation pane.



The Method Development navigation pane opens.

When you save a method, the application adds it to the Recent Files list. The Recent Files list displays a list of your most recently saved master method files.



2. Click the method name in the Recent Files list.

The General page for the selected method opens in the Method View. For a detailed description of all the features on the General page, see "General page" on page 111.

Editing the General Page

The General page defines basic information about the master method. For a detailed description of all the features on the General page, see "General page" on page 111.

Use the following procedures:

- To specify general information for a master method
- To edit the instrument method parameters
- To set automated background subtraction options
- To specify a chromatogram reference sample
- To specify mass tolerance

* To specify general information for a master method

1. In the Lab Name box, type the name to be displayed on the top of each printed, saved, or exported report.

The default name is Default Laboratory.

- 2. In the Assay Type box, type the assay type to be targeted by the method.
- 3. From the Injection Volume box, select the injection volume (in μL) to be used for sample injection.

Use either the up/down arrows to change the volume in increments/decrements of 1 μL or the keyboard to enter non-integer injection volumes.

IMPORTANT The TraceFinder application uses this injection volume in the master method, not the injection volume from the instrument method.

4. From the Ion Range Calc Method list, select a method for calculating the ion ratio range windows.

When you select Level, the TraceFinder application displays a Use Level list where you can choose a calibration level. To define the calibration levels on the Compounds page, see "Editing the Compounds Page" on page 114.

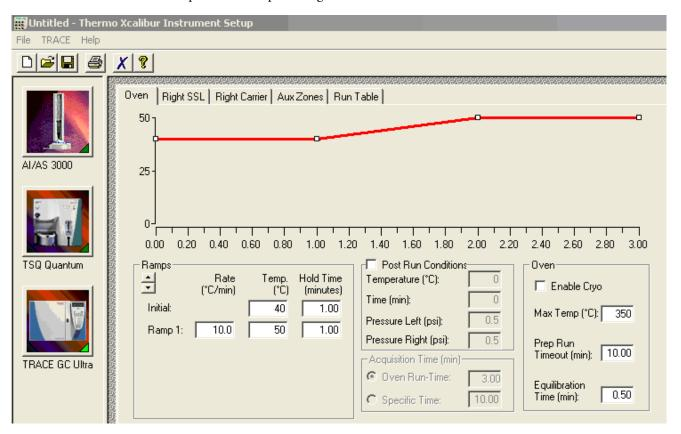


5. From the Qualitative Peak Processing Template box, select a template for performing peak detection on quantitative samples following target compound analysis.

* To edit the instrument method parameters

- From the Instrument Method list on the General page, select an instrument method.
 For a detailed description of all the features on the General page, see "General page" on page 111.
- 2. To edit the instrument method for this master method, click **Edit**.

The Thermo Xcalibur Instrument Setup dialog box opens. This example instrument setup shows multiple configured instruments.

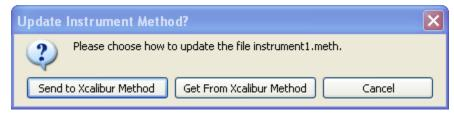


- 3. Edit the values on the instrument page for your instrument.
- 4. From the main menu on the Thermo Xcalibur Instrument Setup dialog box, choose **File > Save** and then choose **File > Exit**.

The TraceFinder application returns you to the General page.

5. To update any changes that were made to the instrument method after you created this master method, click **Update**.

The Update Instrument Method? dialog box opens.



- 6. Choose one of the following options:
 - **Send to Xcalibur Method**: Overwrites the Xcalibur method with the current instrument method.
 - **Get From Xcalibur Method**: Overwrites the current instrument method with the Xcalibur method.
 - Cancel: Make no changes to the instrument method in the current master method.

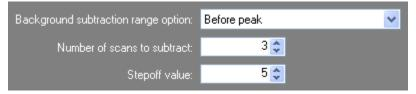
To set automated background subtraction options

- 1. In the Background Subtraction Range Option list, select how you want the subtraction range determined from the following options:
 - **Before Peak**: Averages and subtracts a specified number of scans before the apex of the peak.
 - After Peak: Subtracts a specified number of scans following the apex of the peak.
 - **Both Sides of Peak**: Subtracts a specified number of scans from each side of the apex of the peak.
- 2. In the **Number of Scans to Subtract** box, enter a number.

This is the number of scans that the TraceFinder application subtracts from the background after averaging. If you specified to subtract scans from both sides of the peak, the application subtracts this number of scans from **each** side of the peak.

3. In the Stepoff Value box, enter a number.

The TraceFinder application uses this offset value to average and subtract scans that are not adjacent to the apex of the peak. For example:

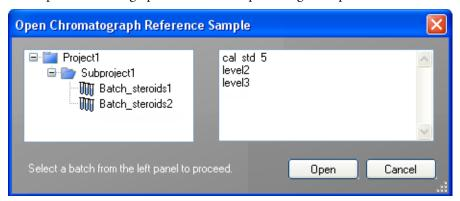


If you specified to subtract 3 scans before the peak and the stepoff value is 5, the TraceFinder application ignores the first 5 scans to the left of the peak and applies the averaging and subtraction to the 6th, 7th, and 8th scans to the left of the peak.

❖ To specify a chromatogram reference sample

- 1. In the Set Chromatogram Reference Sample list, select External.
- 2. Click Select.

The Open Chromatograph Reference Sample dialog box opens.



Note If you are creating a new method, you will not see any reference samples here. You must first create and save a batch using this 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 TraceFinder application displays only batches that were created using the current master method.

6. Select a sample from the list of processed samples.

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

7. Click Open.

The selected sample is displayed as the chromatogram reference sample in the Master Method View.

Tip To clear the reference sample from the master method, select **None** in the Set Chromatogram Reference Sample list.

4 Using the Method Development Mode

Working with Master Methods

❖ To specify mass tolerance

- 1. Select the units of measure you want to use.
- 2. Specify the number of millimass units or parts per million to use as the m/z ± tolerance value.

The application applies this mass tolerance to the extracted chromatograms.

Figure 22. General page

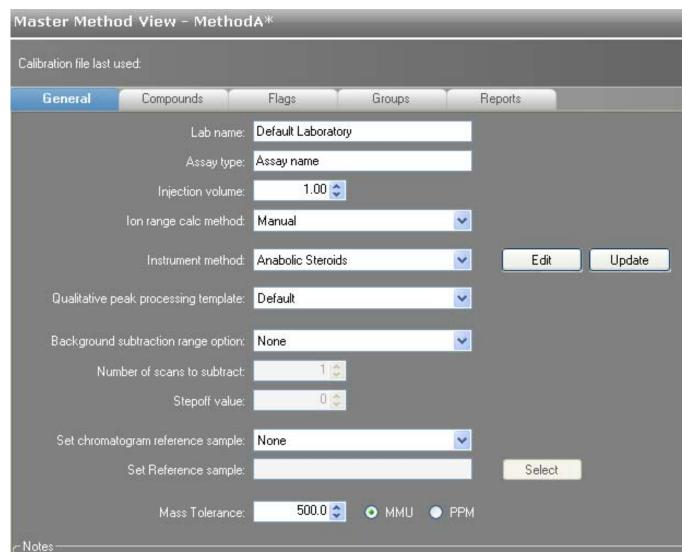


Table 21. General page parameters (Sheet 1 of 3)

| Parameter | Description |
|------------|--|
| Lab name | The laboratory name to be displayed on the top of each printed, saved, or exported report. Default: Default Laboratory To specify this default laboratory name, see "Specifying Configuration Defaults" on page 68. |
| Assay type | The name for the analysis type to be targeted by the method. The assay type associates the method with the analysis of a compound or specific class of compounds (for example, you might use an assay type of PAH for the analysis of Polynuclear Aromatic Hydrocarbons). The TraceFinder application uses this assay type in the method template. |

4 Using the Method Development Mode Working with Master Methods

Table 21. General page parameters (Sheet 2 of 3)

| Parameter | Description |
|---|--|
| Injection volume | The system uses the injection volume (in μL) for sample injection. For a more detailed explanation, refer to the documentation for the autosampler. |
| | The injection volume in the master method overrides the injection volume in the instrument method. |
| | The injection volume in the batch overrides the injection volume in the master method. |
| | Range: 0.1 through 5000 µL |
| Ion range calc method | The TraceFinder application uses the selected ion range calc method to calculate the ion ratio range windows: Manual (default), Average, Level, or Weighted average. When you select Level, an additional list is displayed where you can select a calibration level amount. To define these calibration levels on the Compounds page, see "Editing the Compounds Page" on page 114. |
| Instrument method | Instrument method used for acquiring samples. |
| Edit | Opens the Thermo Xcalibur Instrument Setup dialog box where you can edit the instrument method. |
| Update | Choose one of the following: Send to Xcalibur Method: Overwrites the Xcalibur method with the current instrument method. Get From Xcalibur Method: Overwrites the current instrument method with the Xcalibur method. |
| Qualitative peak processing template | The TraceFinder application uses the qualitative peak processing template to perform peak detection on quantitative samples following compound analysis. |
| Background subtraction range option | Valid values: None, Before Peak, After Peak, Both Sides of Peak Default: None |
| Number of scans to subtract | Valid values: Even numbered integers Default: 0 |
| Stepoff value | Offset from the selected peak to the first subtracted peak. |
| Set chromatogram reference sample | Valid values: None, External Default: None |
| Set Reference sample | This parameter is enabled only when Set Chromatogram Reference Sample is set to External. Click the Select button to choose a reference sample from the project folders. |

Table 21. General page parameters (Sheet 3 of 3)

| Parameter | Description |
|----------------|---|
| Mass Tolerance | Upper limit of MMU or PPM. Default: 500 Range: 0.1 through 50 000 |
| | (Default) MMU (millimass units): MMU is a static calculation to the extracted mass. PPM (parts per million): PPM is a variable calculation dependent on the actual mass. The smaller the mass, the narrower the tolerance range. The larger the mass, the wider the tolerance range. |
| Notes | Notes you add about the method. |

Editing the Compounds Page

Use the Compounds page to set all parameters for the identification, detection, and quantification for the target compound list.

From the Compounds page of the Master Method View, you can access the following pages:

- Acquisition List
- Identification
- Detection
- Calibration
- Calibration Levels
- QC Levels
- Real Time Viewer
- Using the Shortcut Menu Commands

Acquisition List

The Acquisition List page displays all compounds defined for the current method. From this page, you can add or delete compounds from the method. For a detailed description of all the features on the Acquisition List page, see "Acquisition List page" on page 116.

The Acquisition List page is displayed only when you select the Enable Compound Datastore option on the Compound Datastore page in the Application Configuration view. For a detailed description of all the features on the Application Configuration view., see "Application Configuration" on page 56.

Use the following procedures:

- To filter the compound list
- To delete a compound from the list
- To add a compound to the list

❖ To filter the compound list

1. To display a filtered list of compounds, click the funnel icon, , in the column header.

The application displays a list of filterable criteria. In all lists, you can choose to filter by All, Blanks, NonBlanks, or by custom filter criteria. Other filter criteria are specific to the individual columns.

2. To create a custom filter based on compound values in a specific column, choose **Custom** from the column list.

For detailed instructions about creating a custom filter, see Appendix C, "Using Filter Criteria."

❖ To delete a compound from the list

- 1. Select the compound to remove from the list.
- Click the Remove Compound icon, or right-click and choose Remove Compound from the shortcut menu.

A confirmation dialog box opens, listing the compound to be removed.

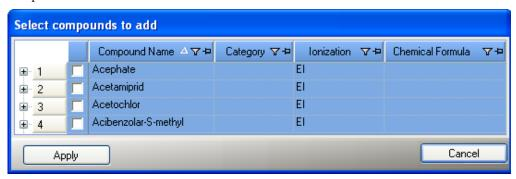
3. To confirm the deletion, click **Yes**.

The selected compound is removed from the acquisition list, which has no effect on the compound datastore.

To add a compound to the list

1. Click the **Add Compound** icon, , or right-click and choose **Add Compound** from the shortcut menu.

The Select Compounds to Add dialog box opens, listing all the compounds defined in the compound datastore.



- 2. Select the check box for each of the compounds you want to add to the method.
- 3. To quickly select multiple compounds, right-click and use any of the commands on the shortcut menu.

Table 22. Select Compounds to Add shortcut menu

| Command | Description |
|--------------|--|
| Select All | Selects all compounds in the compound datastore. |
| Deselect All | Clears all compounds in the compound datastore. |
| Copy Down | Copies the state (checked or cleared) of the currently selected compound to all compounds below it in the compound list. |

4. Click Apply.

The TraceFinder application adds the compounds to the Acquisition List and Compounds pages of the Master Method View.

4 Using the Method Development Mode

Working with Master Methods

Figure 23. Acquisition List page

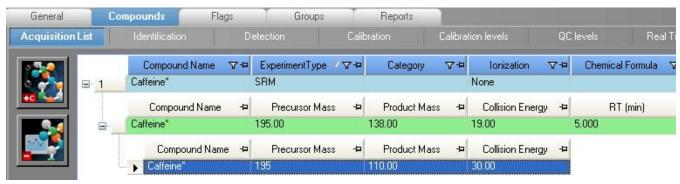


Table 23. Acquisition List page parameters (Sheet 1 of 3)

| Parameter | Description | |
|----------------------|---|--|
| Function Icons | | |
| *** | Opens the Select Compounds to Add dialog box that lists all the compounds defined in the compound datastore. | |
| | Deletes the selected compound. The icon is unavailable when no row is selected. If you used the filters to display a subset of compounds, the selected compound might not be visible on the Acquisition List page. | |
| Compound parameter | | |
| Compound Name | Alphanumeric name assigned to the compound. | |
| Experiment Type | Experiment type: SRM, XIC, or SIM. | |
| Category | (Optional) Alphanumeric identifier. | |
| Ionization | (Optional) Alphanumeric identifier. Valid values: ESI, APCI, EI, CI, APPI | |
| Chemical Formula | (Optional) Alphanumeric chemical identifier. | |
| Quan peak parameters | | |
| Precursor Mass | The mass-to-charge ratio of a precursor ion. The location of the center of a target precursor-ion peak in mass-to-charge ratio (m/z) units. Available only for SRM experiments. Range: 10.000 to 2999.999 | |
| Product Mass | The mass-to-charge ratio of the quan ion. The location of the center of a target quan-ion peak in mass-to-charge ratio (m/z) units. Available only for SRM experiments. Range: 10.000 to 2999.999 | |

Table 23. Acquisition List page parameters (Sheet 2 of 3)

| Parameter | Description | | |
|----------------------|---|--|--|
| Mass | The mass-to-charge ratio of the quan ion. The location of the center of a target quan-ion peak in mass-to-charge ratio (m/z) units. Available only for XIC and SIM experiments. Range: 10.000 to 2999.999 | | |
| Collision Energy | The energy used when ions collide with the collision gas. Range: -250 to 250 | | |
| Lens | (Optional) Range: –400 to 400 | | |
| Polarity | + (positive) or – (negative) | | |
| RT (min) | Retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. | | |
| | The TraceFinder application uses the RT and Window values to determine the start and stop time for the acquisition. Range: 0.00 to 999.00 Start time = RT - (Window/2) Stop time = RT + (Window/2) Start and stop range: 0.00 to 999.00 | | |
| Window (sec) | Acquisition window. The TraceFinder application uses the RT and Window values to determine the start and stop time for the acquisition. Range: 0.00 to 499.50 Start time = RT - (Window/2) Stop time = RT + (Window/2) Start and stop range: 0.00 to 999.00 | | |
| Energy Ramp | Available only for SRM experiments. Range: 0.00 to 200.00 | | |
| Confirming ion param | neters | | |
| Precursor Mass | The mass-to-charge ratio of a precursor ion. The location of the center of a target precursor-ion peak in mass-to-charge ratio (<i>m/z</i>) units. Available as a read-only field for SRM experiments only. Range: 10.000 to 2999.999 | | |
| Product Mass | The mass-to-charge ratio of the quan ion. The location of the center of a target quan-ion peak in mass-to-charge ratio (m/z) units. Available only for SRM experiments. Range: 10.000 to 2999.999 | | |

4 Using the Method Development Mode Working with Master Methods

Table 23. Acquisition List page parameters (Sheet 3 of 3)

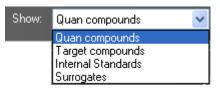
| Parameter | Description |
|------------------|---|
| Mass | The mass-to-charge ratio of the quan ion. The location of the center of a target quan-ion peak in mass-to-charge ratio (m/z) units. Available only for XIC and SIM experiments. Range: 10.000 to 2999.999 |
| Collision Energy | The energy used when ions collide with the collision gas. Available only for SRM experiments. Range: -250.00 to 250.00 |

Identification

The Identification page lists the compounds that are targeted for analysis, reporting, and other compound-specific values.

❖ To filter the displayed compounds

From the Show list, select the type of compounds you want to display in the compounds list.



| Compound type | Description |
|--------------------|---|
| Quan compounds | Displays only quan compounds, such as target compounds, internal standards, and surrogates. |
| Target compounds | Displays only target compounds. |
| Internal Standards | Displays only internal standard compounds. |
| Surrogates | Displays only surrogate compounds. |

Figure 24. Identification page

| General | | Compounds | Flags G | Groups | Repo | orts | |
|-------------|------|---------------------|-------------------|--------|------------|---------------------|--------------------|
| Acquisition | List | Identification | Detection | С | alibration | Calibration levels | s QC levels |
| | RT | Compound | Compound type | Active | CAS No | Use as RT Reference | Reference compound |
| 1 | 4.38 | 1,3-Dichlorobenzene | Target Compound | ~ | | | |
| 2 | 4.45 | 1,4-Dichlorobenzene | Internal Standard | ~ | | V | |

Table 24. Identification page parameters (Sheet 1 of 2)

| Parameter | Description | | |
|---|---|--|--|
| RT Retention time. The time after injection when the compound elutes. The total t compound is retained on the column. | | | |
| | The TraceFinder application uses the RT and Window values to determine the start and stop time for the acquisition. Range: 0.00 to 999.00 Start time = RT - (Window/2) Stop time = RT + (Window/2) Start and stop range: 0.00 to 999.00 | | |
| Compound | A list of compounds that have been identified. To customize the compound names, click the cell and type a new name. To display a filtered list of compounds, use the Show list. | | |

4 Using the Method Development Mode Working with Master Methods

Table 24. Identification page parameters (Sheet 2 of 2)

| Parameter | Description | | | |
|---------------------|--|--|--|--|
| Compound type | Compound types are Quan Compound, Target Compound, Internal Standard, and Surrogate. | | | |
| | The TraceFinder application uses target compounds, internal standards, and surrogates in quantitative analysis. | | | |
| Active | Identifies each compound to be included in data review and reporting. By default, all added compounds are set to active. This active or inactive setting populates to the Batch View and Data Review view in the Data Review mode. | | | |
| CAS No | The Chemical Abstract Service (CAS) number that the TraceFinder application matched with each compound. To change or add a number, click the CAS No cell and enter a new number. | | | |
| Use as RT Reference | When performing peak detection with retention time standards, the TraceFinder application first identifies those compounds identified as retention time standards and then uses their observed retention times to adjust any associated target compound. | | | |
| Reference compound | Reference compound to be used for retention time adjustment for a compound. This list includes all compounds that are selected in the Use as RT Reference column. | | | |

Detection

Use the Detection page to customize peak detection and integration for any ions that define peaks and compounds.

From the Detection page, you can access the following pages:

- Times
- Signal
- Detect
- Spectrum
- Ratios

On the Detection page (see "Detection page" on page 129), you can configure how characteristic ions for targeted compounds are detected and integrated. You can also edit the list of characteristic ions for a specific compound. Refining these parameters in the master method for each compound and its ions can reduce the degree of manual integration that would otherwise be required.

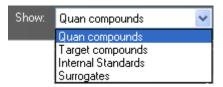
You can change the parameters used to identify a quantification peak, mass range, or confirming ion. The TraceFinder application automatically uses the first match it finds as the compound name, the base peak of the mass spectrum as the quan peak, and the second and third largest ions as the confirming ions.

Use the following procedures:

- To filter the displayed compounds
- To add compounds to the method
- To change the compound reference spectrum
- To replace a quan mass
- To add a mass to the existing quan mass ranges
- To add a quan peak
- To add a mass as a new compound
- To replace a confirming ion
- To add a mass as a new confirming ion
- To save the new method

❖ To filter the displayed compounds

From the Show list, select the type of compounds you want to display in the compounds list.

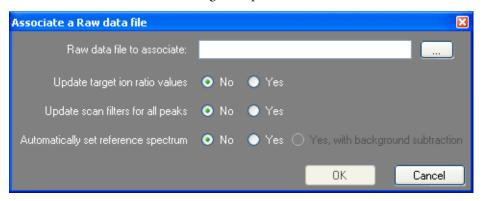


| Compound type | Description |
|--------------------|--|
| Quan compounds | Displays all quan compounds, such as target compounds, internal standards, and surrogates. |
| Target compounds | Displays only target compounds. |
| Internal Standards | Displays only internal standard compounds. |
| Surrogates | Displays only surrogate compounds. |

To add compounds to the method

1. From the main menu, choose Master Method > Associate a Raw Data File.

The Associate a Raw Data File dialog box opens.



- 2. Browse to a raw data file to associate with the method and open the file.
- 3. To update the target ion ratio values when you associate this raw data file, select the **Yes** option.
- 4. To update the scan filters when you associate this raw data file, select the **Yes** option.
- To set a reference spectrum, do one of the following:Select the **Yes** option.

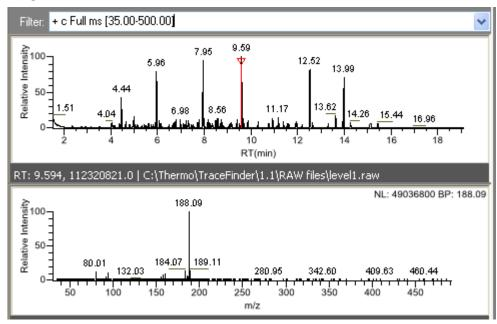
-Or-

Select the Yes, with Background Subtraction option.

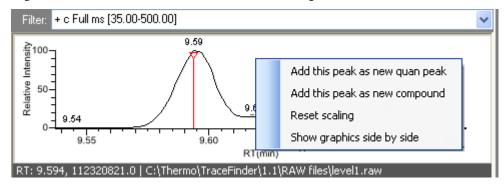
This feature is available only when you have set background subtraction values on the General page of the Master Method View. See "Editing the General Page" on page 106.

6. Click OK.

The TraceFinder application displays the chromatographic and spectrum data for the compounds in the selected raw data file.



- 7. Select a filter from the Filter list.
- 8. Click to select the peak in the chromatogram that represents the compound you want to add to the method.
- 9. Right-click and choose Add This Peak as New Compound from the shortcut menu.



The TraceFinder application performs a library search for the selected compound. The application uses the first match it finds as the compound name, the base peak of the mass spectrum as the quan peak, and the second and third largest ions as the confirming ions.

If the name of the first match is already in the library, the Add New Compound dialog box opens.

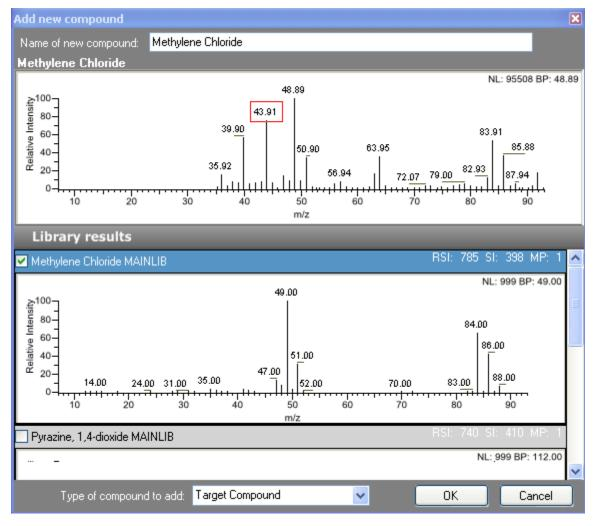
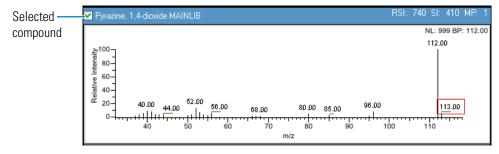


Figure 25. Add New Compound dialog box

10. (Optional) To use a compound other than the compound already in the library, scroll to the spectrum for that compound and select the compound name in the title bar of the spectrum pane.



- 11. In the Type of Compound to Add list, select a compound type.
- 12. Click **OK**.
- 13. Repeat this procedure for each compound you want to add to the method.

To change the compound reference spectrum

1. In the raw data file chromatogram pane, click a peak.

The TraceFinder application displays the spectrum for the selected peak in the spectrum pane.

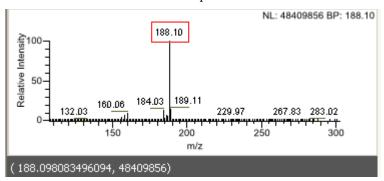
2. In the raw data file spectrum pane, right-click and choose **Use This Spectrum for Compound Reference Spectrum** from the shortcut menu.

The TraceFinder application replaces the spectrum on the Spectrum page of the quan peak pane with this spectrum.

To replace a quan mass

- 1. Click the pane for the quan mass that you want to replace.
- 2. In the raw data file spectrum pane, hold the cursor over a peak.

The red box indicates the selected peak.



- 3. Right-click and choose **Set This Mass as Quan Mass** from the shortcut menu.
- 4. Choose either Don't Update Ion Ratios or Update Ion Ratios Using This Spectrum.

You can see the updated ion ratios on the Ratios page for the confirming ions. See "Ratios" on page 149.

❖ To add a mass to the existing quan mass ranges

1. In the raw data file spectrum pane, hold the cursor over a peak.

The red box indicates the selected peak.

- 2. Right-click and choose **Add This Mass to Existing Quan Mass Ranges** from the shortcut menu.
- 3. Choose either Don't Update Ion Ratios or Update Ion Ratios Using This Spectrum.

The TraceFinder application adds the selected mass to the existing quan mass ranges to increase the signal.

If you chose to update the ion ratios, you can see the updated ion ratios on the Ratios page for the confirming ions. See "Ratios" on page 149.

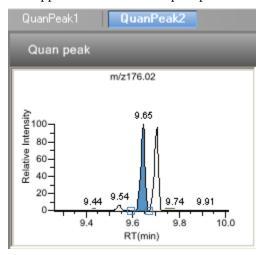
To add a quan peak

1. In the raw data file spectrum pane, hold the cursor over a peak.

The red box indicates the selected peak.

2. Right-click and choose Add This Mass as New Quan Peak from the shortcut menu.

The application adds a new quan peak to the compound.



You can use the shortcut menu on the spectrum pane for this new quan peak to perform any of the tasks you could perform on the original quan peak.

To add a mass as a new compound

1. In the raw data file spectrum pane, hold the cursor over a peak.

The red box indicates the selected peak.

2. Right-click and choose Add This Mass as New Compound from the shortcut menu.

The TraceFinder application performs a library search for the selected compound. The application uses the first match it finds as the compound name, the base peak of the mass spectrum as the quan peak, and the second and third largest ions as the confirming ions.

If the name of the first match is already in the library, the Add New Compound dialog box opens with the matching compound selected.

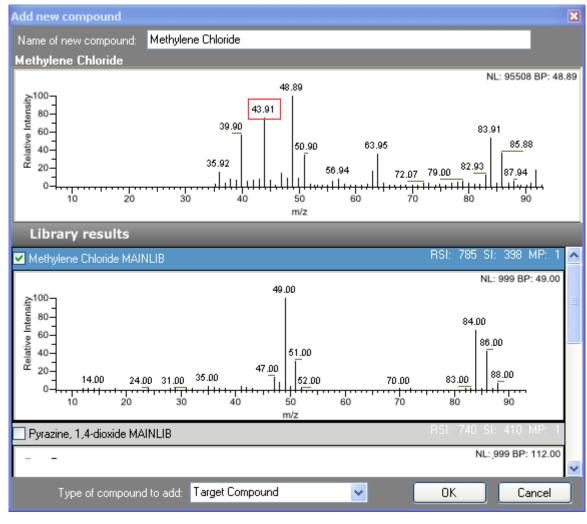
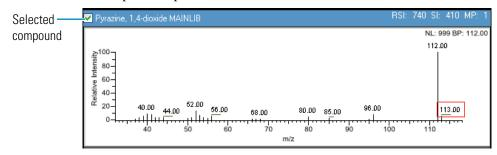


Figure 26. Add New Compound dialog box

- 3. (Optional) In the Add New Compound dialog box, make any of the following changes:
 - a. Change the name for the compound in the Name of New Compound box.
 - b. Use a compound other than the compound chosen by the TraceFinder application by scrolling to the spectrum for that compound and selecting the compound name in the title bar of the spectrum pane.



- c. In the Type of Compound to Add list, select a compound type.
- 4. Click OK.

4 Using the Method Development Mode

Working with Master Methods

To replace a confirming ion

- 1. Click the pane for the confirming ion that you want to replace.
- 2. In the raw data file spectrum pane, hold the cursor over a peak.

The red box indicates the selected peak.

3. Right-click and choose **Set this Mass as Confirming Ion** from the shortcut menu. The TraceFinder application replaces the confirming ion with the selected mass.

❖ To add a mass as a new confirming ion

- In the raw data file spectrum pane, hold the cursor over a peak.
 The red box indicates the selected peak.
- 2. Right-click and choose **Add this Mass as New Confirming Ion** from the shortcut menu. The TraceFinder application adds the confirming ion to the quan peak.



You can use the shortcut menu on the spectrum pane for this new confirming ion to perform any of the tasks you could perform on the original confirming ions.

❖ To save the new method

1. Choose **File > Save**.

The Save Master Method dialog box opens.

2. Do one of the following:

Type a new name for the master method and click **OK**.

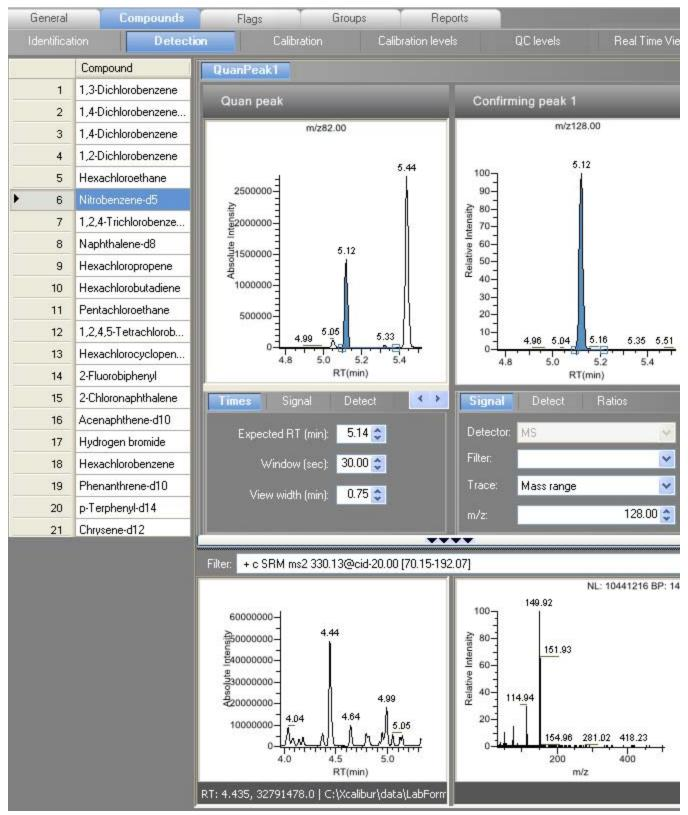
-Or-

Select a method name to overwrite and click **Overwrite**.

The TraceFinder application saves the new method data in the following folder:

...\Thermo\TraceFinder\1.1\Methods

Figure 27. Detection page



4 Using the Method Development Mode Working with Master Methods

 Table 25.
 Detection page parameters

| Parameter | Description |
|--|---|
| Compound list | Lists all compounds in the master method. |
| Quan peak | Displays a chromatogram for the quan peak and its confirming ions. The quan peak and confirming ion panes include additional pages for retention time, signal, detection, spectrum, and ratio parameters. |
| Filter | Displays the filter used for the raw data file. |
| Reference chromatogram and spectra | Displays a reference chromatogram and spectra for the raw data file. |
| Additional pages | |
| Times | Defines the retention time and window for a quan peak. See "Times" on page 131. |
| Signal | Defines the detector and its parameters used to display each chromatogram trace. See "Signal" on page 132. |
| Detect | Defines the peak detection algorithm and its options. See "Detect" on page 135. |
| Spectrum | Defines a reference mass spectrum for a quan peak or compound. See "Spectrum" on page 144. |
| Ratios | Defines the criteria for evaluating, confirming, or qualifying ions. See "Ratios" on page 149. |

Times

Use the Times page to define the retention time and window for a quan peak.

Figure 28. Times page

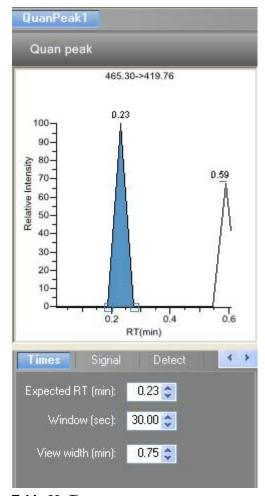


Table 26. Times page parameters

| Parameter | Description |
|-------------------|---|
| Expected RT (min) | Expected retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| Window (sec) | Width of the window (in seconds) to indicate how far around the expected retention time the system looks for a peak apex. |
| View width (min) | Viewable size of the ion chromatogram display. Changing the view width does not affect the process of peak detection; the TraceFinder application uses it only for graphical display. |

Signal

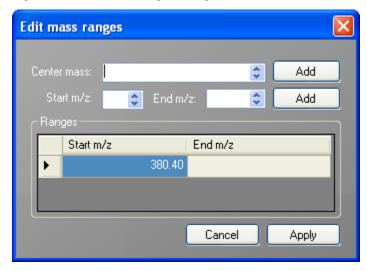
Use the Signal page to define the detector and its parameters as you display each chromatogram trace. For a detailed description of all the features on the Signal page, see "Signal page parameters" on page 133.

To specify ranges of ions for detection and integration

1. In the Ranges area, click Edit.

The Edit Mass Ranges dialog box opens where you can enter rows using a center of mass or a range.

Figure 29. Edit Mass Ranges dialog box



2. Enter a value in the Center Mass box and click Add.

The application adds a row with this value to the ranges grid. Center mass values are listed in the Start m/z column. The application uses a range of one amu centered on this value.

3. Enter values in the Start m/z and End m/z columns and click **Add**.

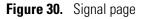
The application adds a row with these start and end values.

4. Add as many ranges as you want.

When you process a batch with this method, the application sums the multiple ions specified by these ranges.

5. Click Apply.

When you are editing a master method, the application applies the entries to the list of ranges.



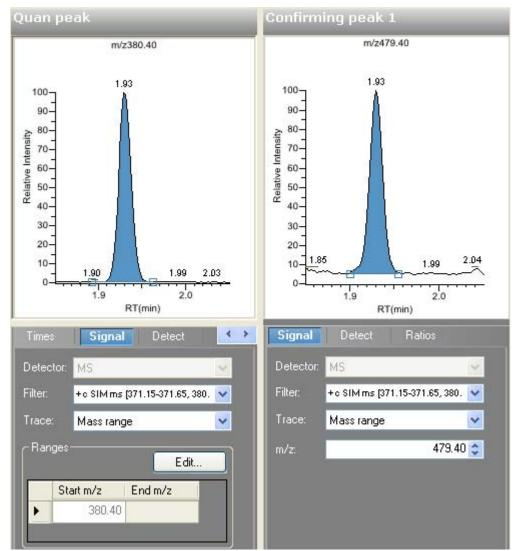


Table 27. Signal page parameters (Sheet 1 of 2)

| Parameter | Description |
|-----------|---|
| Detector | MS: Mass spectrometers. Analog: Supplemental detectors (for example, FID, ECD). AD card: If you have a detector not under data system control, you can capture the analog signal and convert it to digital using an interface box (for example, SS420X) for storage in the raw data file. |
| Filter | Represents a particular data acquisition channel. For example, the filter option + c Full ms [35.00-500.00] represents a positive ion centroid signal acquired in single-stage, full-scan mode from m/z 35 to 500. |
| Trace | Represents a specific range of the data. In conjunction with the filter, the TraceFinder application uses the trace to identify the characteristic ions for a compound. The options are: Mass Range, TIC, or Base Peak. |

4 Using the Method Development Mode Working with Master Methods

Table 27. Signal page parameters (Sheet 2 of 2)

| Parameter | Description | | | | |
|----------------------|---|--|--|--|--|
| Ranges | | | | | |
| Edit | Opens the Edit Mass Ranges dialog box where you can specify a range of ions for detection and integration. See "Edit Mass Ranges dialog box" on page 132. | | | | |
| Start m/z End m/z | Specifies ranges of ions for detection and integration. The application sums the multiple ions specified by these ranges. | | | | |
| | Ranges specified by a center mass value are listed as a single value in the Start m/z column. The application uses a range of one amu centered on this value. | | | | |

Detect

Use the Detect page to define the peak detection algorithm (sensitivity) and its options and to determine the area under a curve. There are three sensitivity modes: Genesis, ICIS, and Avalon. On this page, you can specify how you want each mode to run.

For a detailed description of all the features on the Detect page for Genesis sensitivity, see "Detect page parameters for Genesis" on page 137.

For a detailed description of all the features on the Detect page for ICIS sensitivity, see "Detect page parameters for ICIS" on page 140.

For a detailed description of all the features on the Detect page for Avalon sensitivity, see "Detect page parameters for Avalon" on page 142.

❖ To apply current peak detection settings to all peaks

- 1. Edit the parameters for either the Quan peak or a Confirming peak.
- 2. Right-click the Detect page and choose **Apply to All Peaks in Method** from the shortcut menu.

The application updates all compounds in the method with the current settings on the Detect page. These updates apply to both quan and confirming ions.

To apply current peak detection settings to all peaks that use the same detection mode

- 1. Edit the parameters for either the Quan peak or a Confirming peak.
- 2. Right-click the Detect page and choose Apply to All Peaks with Like Sensitivity Setting from the shortcut menu.

The application uses the current settings on the Detect page to update all compounds in the method that use the same sensitivity mode (Genesis, ICIS, or Avalon). These updates apply to both quan and confirming ions that use the same sensitivity mode.

4 Using the Method Development Mode

Working with Master Methods

Figure 31. Detect page for Genesis

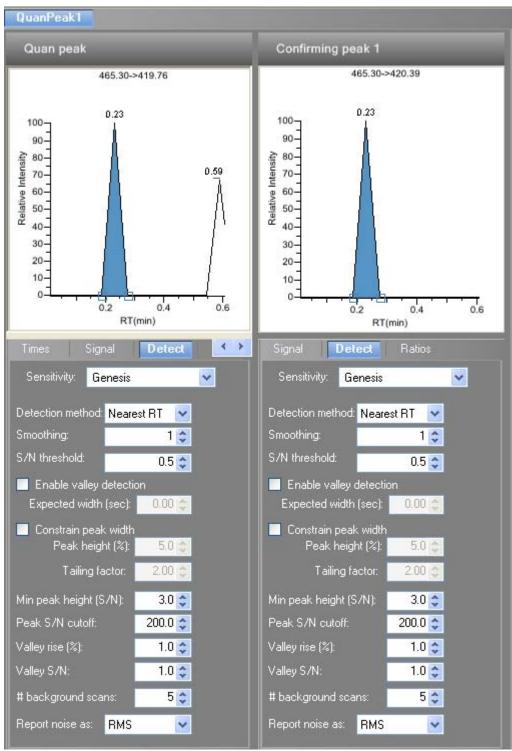


Table 28. Detect page parameters for Genesis (Sheet 1 of 2)

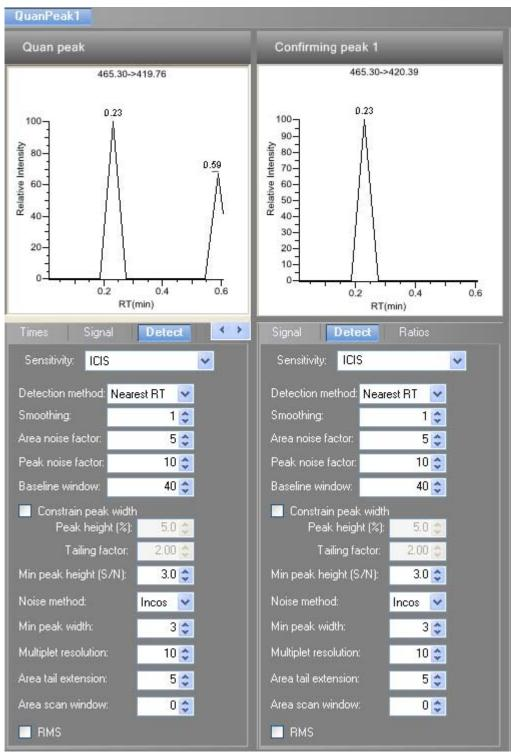
| Parameter | Description | | | |
|---|---|--|--|--|
| Sensitivity Specifies the Genesis peak detection algorithm. | | | | |
| Detection method | Highest peak: Uses the highest peak in the chromatogram for component identification. | | | |
| | Nearest RT: Uses the peak with the nearest retention time in the chromatogram for component identification. | | | |
| Smoothing | Determines the degree of data smoothing to be performed on the active component peak prior to peak detection and integration. | | | |
| | Range: Any odd integer from 1 through 15 points Default: 1 | | | |
| S/N threshold | Current signal-to-noise threshold for peak integration. Peaks with signal-to-noise values less than this value are not integrated. Peaks with signal-to-noise values greater than this value are integrated. | | | |
| | Range: 0.0 to 999.0 | | | |
| Enable valley detection | Uses the valley detection approximation method to detect unresolved peaks. This method drops a vertical line from the apex of the valley between unresolved peaks to the baseline. The intersection of the vertical line and the baseline defines the end of the first peak and the beginning of the second peak. | | | |
| Expected width (sec) | The expected peak width parameter (in seconds). This parameter controls the minimum width that a peak is expected to have if valley detection is enabled. | | | |
| | With valley detection enabled, any valley points nearer than the <i>expected width</i> /2 to the top of the peak are ignored. If a valley point is found outside the expected peak width, the TraceFinder application terminates the peak at that point. The application always terminates a peak when the signal reaches the baseline, independent of the value set for the expected peak width. | | | |
| | Range: 0.0 to 999.0 | | | |
| Constrain peak width | Constrains the peak width of a component during peak integration of a chromatogram. You can then set values that control when peak integration is turned on and off by specifying a peak height threshold and a tailing factor. Selecting the Constrain Peak Width check box enables the Peak Height (%) and Tailing Factor options. | | | |
| Peak height (%) | A signal must be above the baseline percentage of the total peak height (100%) before integration is turned on or off. This text box is active only when you select the Constrain Peak Width check box. | | | |
| | | | | |

4 Using the Method Development Mode Working with Master Methods

Table 28. Detect page parameters for Genesis (Sheet 2 of 2)

| Parameter | Description | | | | |
|---|---|--|--|--|--|
| Tailing factor | A factor that controls how the TraceFinder application integrates the tail of a peak. This factor is the maximum ratio of the trailing edge to the leading side of a constrained peak. This text box is active only when you select the Constrain the Peak Width check box. | | | | |
| | Range: 0.5 through 9.0 | | | | |
| Min peak height (S/N) | For the valley detection approximation method to use the Nearest RT Peak Identification criteria, this peak signal-to-noise value must be equaled or exceeded. For component identification purposes, the TraceFinder application ignores all chromatogram peaks that have signal-to-noise values that are less than the S/N Threshold value. | | | | |
| | Range: 0.0 (all peaks) through 999.0 | | | | |
| Peak S/N cutoff | The peak edge is set to values below this signal-to-noise ratio. | | | | |
| | This test identifies an edge of a peak when the baseline adjusted height of the edge is less than the ratio of the baseline adjusted apex height and the peak S/N cutoff ratio. | | | | |
| | When the S/N at the apex is 500 and the peak S/N cutoff value is 200, the TraceFinder application defines the right and left edges of the peak when the S/N reaches a value less than 200. | | | | |
| | Range: 50.0 to 10000.0 | | | | |
| Valley rise (%) | The peak trace can rise above the baseline by this percentage after passing through a minimum (before or after the peak). | | | | |
| | This method drops a vertical line from the apex of the valley between unresolved peaks to the baseline. The intersection of the vertical line and the baseline defines the end of the first peak and the beginning of the second peak. | | | | |
| | When the trace exceeds rise percentage, the TraceFinder application applies valley detection peak integration criteria. | | | | |
| | The TraceFinder application applies this test to both the left and right edges of the peak. | | | | |
| | The rise percentage criteria is useful for integrating peaks with long tails. | | | | |
| | Range: 0.1 to 500.0 | | | | |
| Valley S/N | Specifies a value to evaluate the valley bottom. Using this parameter ensures that the surrounding measurements are higher. | | | | |
| | Range: 1.0 to 100.0 Default: 2.0 | | | | |
| # background scans | Number of background scans performed by the TraceFinder application. | | | | |
| Report noise as Determines if the noise used in calculating S/N values is calculated using an RM calculation or a peak-to-peak resolution threshold. Options are RMS or Peak to I | | | | | |

Figure 32. Detect page for ICIS



4 Using the Method Development Mode Working with Master Methods

Table 29. Detect page parameters for ICIS (Sheet 1 of 2)

| Description | | | |
|--|--|--|--|
| Specifies the ICIS peak detection algorithm. | | | |
| Highest peak: Uses the highest peak in the chromatogram for component identification. | | | |
| Nearest RT: Uses the peak with the nearest retention time in the chromatogram for component identification. | | | |
| Determines the degree of data smoothing to be performed on the active component peak prior to peak detection and integration. | | | |
| Range: Any odd integer from 1 through 15 points Default: 1 | | | |
| The noise level multiplier used to determine the peak edge after the location of the possible peak. | | | |
| Range: 1 through 500 Default: 5 | | | |
| The noise level multiplier used to determine the potential peak signal threshold. | | | |
| Range: 1 through 1000 Default: 10 | | | |
| The TraceFinder application looks for a local minima over this number of scans. | | | |
| Range: 1 through 500 Default: 40 | | | |
| Constrains the peak width of a component during peak integration of a chromatogram. You can then set values that control when peak integration is turned on and off by specifying a peak height threshold and a tailing factor. Selecting the Constrain Peak Width check box enables the Peak Height (%) and Tailing Factor options. | | | |
| A signal must be above the baseline percentage of the total peak height (100%) before integration is turned on or off. This text box is active only when you select the Constrain Peak Width check box. | | | |
| Range: 0.0 to 100.0% | | | |
| A factor that controls how the TraceFinder application integrates the tail of a peak. This factor is the maximum ratio of the trailing edge to the leading side of a constrained peak. This text box is active only when you select the Constrain the Peak Width check box. | | | |
| Range: 0.5 through 9.0 | | | |
| | | | |

Table 29. Detect page parameters for ICIS (Sheet 2 of 2)

| Parameter | Description | | |
|-----------------------|---|--|--|
| Min peak height (S/N) | For the valley detection approximation method to use the Nearest RT Peak Identification criteria, this peak signal-to-noise value must be equaled or exceeded. For component identification purposes, the TraceFinder application ignores all chromatogram peaks that | | |
| | have signal-to-noise values that are less than the S/N Threshold value. Range: 0.0 (all peaks) through 999.0 | | |
| Noise method | The options are INCOS or Repetitive. | | |
| | INCOS: Uses a single pass algorithm to determine the noise level. | | |
| | Repetitive: Uses a multiple pass algorithm to determine the noise level. In general, this algorithm is more accurate in analyzing the noise than the INCOS Noise algorithm, but the analysis takes longer. | | |
| Min peak width | The minimum number of scans required in a peak. | | |
| | Range: 0 to 100 scans Default: 3 | | |
| Multiplet resolution | The minimum separation in scans between the apexes of two potential peaks. This is a criterion to determine if two peaks are resolved. | | |
| | Range: 1 to 500 scans Default: 10 | | |
| Area tail extension | The number of scans past the peak endpoint to use in averaging the intensity. | | |
| | Range: 0 to 100 scans Default: 5 | | |
| Area scan window | The number of allowable scans on each side of the peak apex. A zero value defines all scans (peak-start to peak-end) to be included in the area integration. | | |
| | Range: 0 to 100 scans Default: 0 | | |
| RMS | Specifies that the TraceFinder application calculate noise as RMS. By default, the application uses Peak To Peak for the noise calculation. RMS is automatically selected if you manually determine the noise region. | | |

4 Using the Method Development Mode

Working with Master Methods

Figure 33. Detect page for Avalon

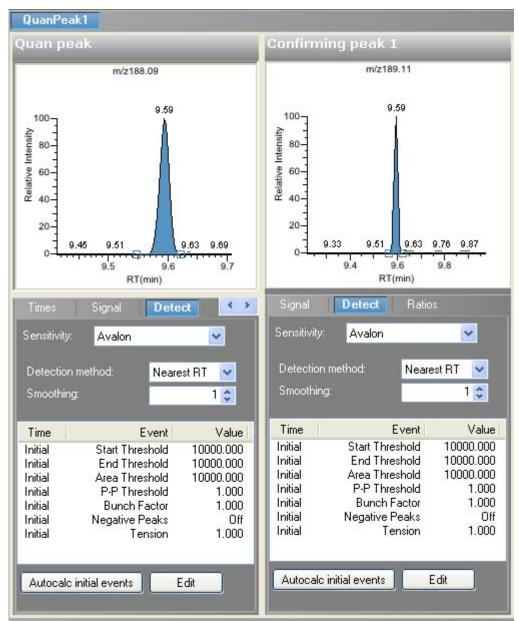


Table 30. Detect page parameters for Avalon (Sheet 1 of 2)

| Parameter | Description | | | | |
|--|---|--|--|--|--|
| Sensitivity Specifies the Avalon peak detection algorithm. | | | | | |
| Detection method | Highest Peak: Uses the highest peak in the chromatogram for component identification. | | | | |
| | Nearest RT: Uses the peak with the nearest retention time in the chromatogram for component identification. | | | | |

Table 30. Detect page parameters for Avalon (Sheet 2 of 2)

| Parameter | Description | | | |
|-------------------------|---|--|--|--|
| Smoothing | Determines the degree of data smoothing to be performed on the active component peak prior to peak detection and integration. | | | |
| | Range: Any odd integer from 1 through 15 points Default: 1 | | | |
| Autocalc Initial Events | Automatically calculates the events in the Event list. | | | |
| Edit | Opens the Avalon Event List dialog box. See "Avalon Event List" on page 80. | | | |

Spectrum

Use the Spectrum page to store a reference mass spectrum for a quan peak or compound.

For a detailed description of all the shortcut menu commands on the Spectrum page, see "Spectrum shortcut menu functions" on page 148.

Use the following procedures:

- To update confirming ion ratios
- To change the quantitation mass used for a quan peak
- To add ions together to get an accumulated signal
- To add a quan peak to an existing compound
- To add one or more confirming ions to an existing compound
- To zoom in on the chromatogram or spectrum displays

To update confirming ion ratios

1. Click a peak in the quan peak chromatogram pane.

The mass spectrum for the peak is displayed in the Spectrum pane.

2. Right-click the Spectrum pane and choose **Update Confirming Ion Ratios with This Spectrum** from the shortcut menu.

❖ To change the quantitation mass used for a quan peak

1. Click a peak in the chromatogram pane.

The mass spectrum for the peak is displayed in the spectrum pane.

2. In the spectrum pane, hold the cursor over the m/z value for an ion.

A red box around the ion's m/z value indicates that the ion is selected.

- 3. Right-click and choose one of the following commands from the shortcut menu:
 - Set This Mass as Quan Mass > Don't Update Ion Ratios
 - Set This Mass as Quan Mass > Update Ion Ratios Using This Reference Spectrum

The following examples show an original quan peak and a quan peak with an updated quan mass.

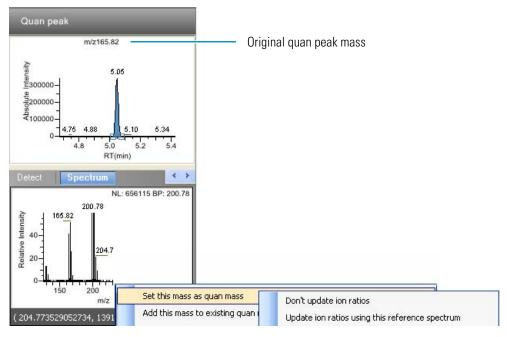


Figure 34. Original quan peak mass example

The TraceFinder application replaces the original quan mass with the selected mass.

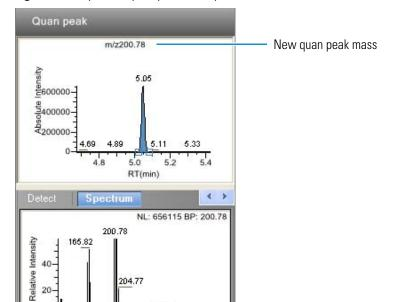


Figure 35. Updated quan peak example

Thermo Scientific TraceFinder User Guide 145

267.02 250

❖ To add ions together to get an accumulated signal

1. Hold the cursor over the m/z value for an ion in the Spectrum pane.

A red box around the ion's *m/z* value indicates that the ion is selected.

2. Right-click and choose **Add This Mass to Existing Quan Mass Range** from the shortcut menu.

You can now update the ion ratios to adjust the confirming ion comparisons to the new summed quan peak signal.

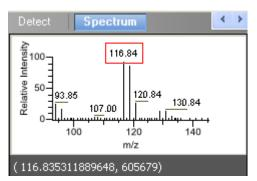
To add a quan peak to an existing compound

1. Click the peak in the Quan Peak chromatogram pane.

The mass spectrum for the peak is displayed in the Spectrum pane.

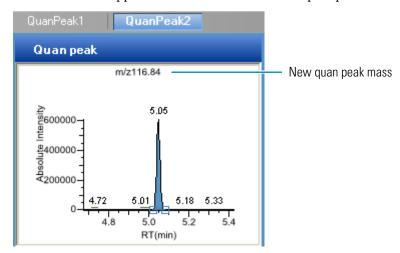
2. In the Spectrum pane, hold the cursor over the m/z value for an ion.

A red box around the ion's m/z value indicates that the ion is selected.



3. Right-click and choose **Set This Mass as New Quan Peak** from the shortcut menu.

The TraceFinder application adds this ion as a new quan peak.



❖ To add one or more confirming ions to an existing compound

1. Click the peak in the chromatogram pane.

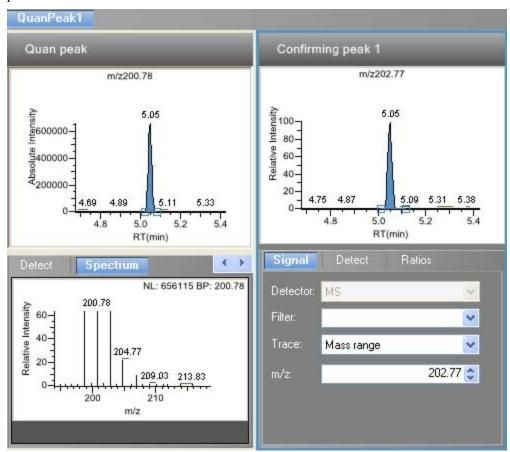
The mass spectrum for the peak is displayed in the Spectrum pane.

2. In the Spectrum pane, hold the cursor over the m/z value for an ion.

A red box around the ion's m/z value indicates that the ion is selected.

3. Right-click and choose to **Add This Mass as New Confirming Ion** from the shortcut menu.

The TraceFinder application adds the selected mass as a confirming peak for this quan peak.



❖ To zoom in on the chromatogram or spectrum displays

1. Drag the cursor to delineate a rectangle.

The display zooms in on the specified rectangle.

2. To return to the original display, right-click and choose **Reset Scaling** from the shortcut menu.

4 Using the Method Development Mode

Working with Master Methods

Figure 36. Spectrum page

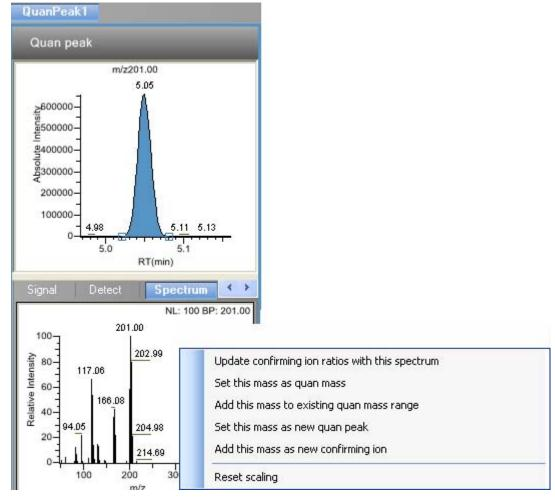


Table 31. Spectrum shortcut menu functions

| Functions | Description | | | |
|---|---|--|--|--|
| Update confirming ion ratios with this spectrum | Updates the confirming ion ratios using the selected peak. | | | |
| Set this mass as quan mass | Adds the quan mass of the selected ion to the quantitation mass used for the quan peak. You can choose to update the ion ratios or not update the ion ratios using this reference spectrum. | | | |
| Add this mass to existing quan mass range | Adds the selected mass to your existing quan mass range. You can choose to update the ion ratios to adjust the confirming ion comparisons to the new summed quan peak signal. | | | |
| Set this mass as new quan peak | Adds a new quan peak to an existing compound. | | | |
| Add this mass as new confirming ion | Adds one or more confirming ions to an existing compound. | | | |
| Reset scaling | Returns the chromatogram or spectrum display to its original size. | | | |

Ratios

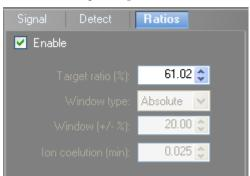
Use the Ratios page to define the criteria for evaluating the confirming or qualifying ions. The TraceFinder application detects compounds that have confirming ion values outside their acceptable window and flags them in the Acquisition mode and on reports.

For a detailed description of all the features on the Ratios page, see "Ratios page parameters" on page 150.

To specify ion ratio criteria

- 1. Select the **Enable** check box to enable the confirming ion.
- 2. In the Target Ratio box, select the theoretical ratio of the confirming ion's response to the quantification ion's response.
- 3. In the Window Type list, select **Absolute** or **Relative** as the calculation approach for determining the acceptable ion ratio range.
- 4. In the Window (+/-%) box, select the acceptable ion ratio range.
- 5. In the Ion Coelution box, select the maximum difference in retention time between a confirming ion peak and the quantification ion peak.

In the following example:



- The target ratio is expected to be 61.02% and the window is Absolute 20%, so the acceptable window for this confirming ion is 41.02 to 81.02%.
- If, instead, the window type is Relative, the plus or minus value is 20% of 61.02% (or 12.20%), so the acceptable window for this confirming ion is 48.82 to 73.22%.

4 Using the Method Development Mode

Working with Master Methods

Figure 37. Ratios page

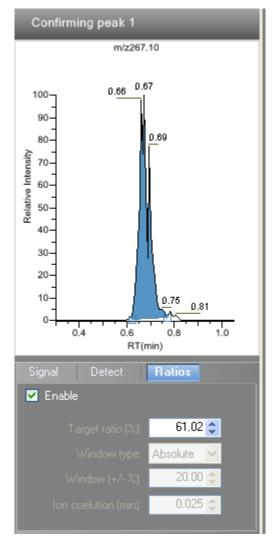


Table 32. Ratios page parameters

| Parameter | Description | | | |
|--|---|--|--|--|
| Enable | Makes the ion ratio criteria available. | | | |
| Target ratio (%) The theoretical ratio of the confirming ion's response to the quantification ion's re | | | | |
| Window type | The absolute or relative calculation approach for determining the acceptable ion ratio range. | | | |
| Window (+/-%) | The acceptable ion ratio range. | | | |
| Ion coelution (min) | The maximum difference in retention time between a confirming ion peak and the quantification ion peak. | | | |

Calibration

Use the Calibration page to set or edit the mathematical model used for preparing the initial calibration evaluation for one or more calibration standards.

Each target compound can have its own initial calibration settings, independent of the other compounds. You can modify the calibration approach on this page or in Acquisition mode when you view the results of an actual calibration batch.

Figure 38. Calibration page

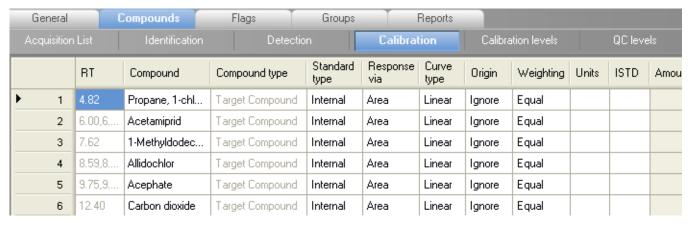


Table 33. Calibration page parameters

| Parameter | Description | | | | |
|---------------|--|--|--|--|--|
| RT | Retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. | | | | |
| Compound | The compound name. | | | | |
| Compound type | Displays the compound type as an internal standard, an external standard, or a target compound. | | | | |
| Standard type | Specifies Internal or External standards. | | | | |
| Response via | The use of area or height. | | | | |
| Curve type | Specifies Linear, Quadratic, or AverageRF curve types. | | | | |
| Origin | The origin treatment as Ignore, Include, or Force. The Origin and Weighting columns are active only when you are using Linear or Quadratic curve types. | | | | |
| Weighting | Specifies the weighting as Equal, 1/X, 1/X ² , 1/Y, or 1/Y ² . | | | | |
| Units | The units to be displayed with the calculated values. | | | | |
| ISTD | The internal standard (ISTD) for a target compound or surrogate if the standard type is set to Internal. If you set the standard type to External, this field is inactive. | | | | |
| Amount | The amount of the internal standard for ISTD compounds. | | | | |

Calibration Levels

On the Calibration levels page for a master method, you can define the standards for calibration. You can edit calibration levels and concentrations for master methods only. The contents of this page are read-only when you are editing a local method.

❖ To specify calibration levels and concentrations

1. Select the compound whose calibration levels and concentrations you want to define.

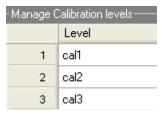


2. In the Manage Calibration Levels area, type a value for the first calibration level.

The TraceFinder application adds a new, empty calibration level row beneath the edited row.



3. Continue adding calibration levels.



When you finish adding calibration levels, you can specify the concentrations for each level for each compound.

- 4. To enter the concentrations to the table, do the following:
 - a. Select the first calibration level table cell.
 - Click the cell again to make it editable.
 - Type a concentration value.

| Acquisition List | | ition List | Identification De | etection | Calibration | Calibration levels | |
|------------------|---|------------|----------------------|----------|-------------|--------------------|------|
| Г | | RT | Compound | cal1 | cal2 | cal3 | cal4 |
| | 1 | 6.00 | Acetamiprid | 10 | | | |
| | 2 | 7.62 | 1-Methyldodecylamine | | | | |

- 5. Repeat Step 4 for all calibration levels associated with the first compound.
- 6. To specify the same concentration values for all compounds, select the value you want to copy, right-click, and choose **Copy Down** from the shortcut menu.

Figure 39. Calibration Levels page

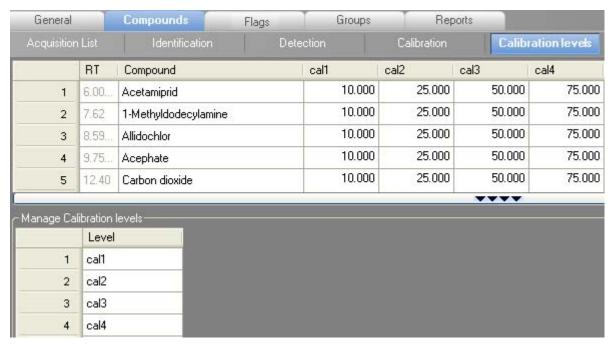


Table 34. Calibration levels page parameters

| Parameter | Description |
|---------------------------|--|
| RT | Retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| Compound | The compound name. |
| cal1-caln | User-defined calibration levels for the compound. |
| Manage Calibration levels | Defines values for each of the calibration level values for the selected compound. |

QC Levels

Use the QC levels page for a master method to define the standards for QC levels. You can edit QC levels for master methods only. The contents of this page are read-only when you are editing a local method. For a detailed description of all the features on the QC Levels page, see "QC levels page parameters" on page 155.

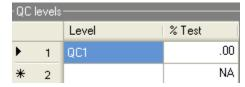
To specify QC levels and concentrations

1. Select the compound whose QC levels, percentage test values, and concentrations you want to define.



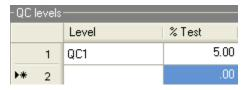
2. In the QC Levels area, type a name for the first QC level.

The TraceFinder application adds a new, empty QC level row beneath the edited row.

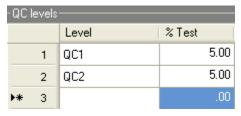


3. Type a value for the % Test.

The % Test is the acceptable difference (as a percentage) between the known amount and the calculated (measured) amount of each QC level.



4. Continue adding QC levels and values for the percentage test.



When you finish adding QC levels, you can specify the concentrations for each level for each compound.

- 5. To enter the concentration values to the table, do the following:
 - a. Select the first QC level table cell.
 - b. Click the cell again to make it editable.
 - c. Type a concentration value.

6. Repeat Step 5 for all QC levels associated with the first compound.

| | Acquisition | List Iden | tification Detection Calibration | Calibration le | vels | QC levels |
|---|-------------|-----------------|----------------------------------|----------------|--------|-----------|
| ľ | | RT | Compound | QC1 | QC2 | QC3 |
| ĺ | 1 | 6.00,6.00,6.00, | Acetamiprid | 10.000 | 15.000 | 25.000 |
| ı | 2 | 7.62 | 1-Methyldodecylamine | | | |

7. To specify the same concentration values for all compounds, select the value you want to copy, right-click, and choose **Copy Down** from the shortcut menu.

Figure 40. QC Levels page

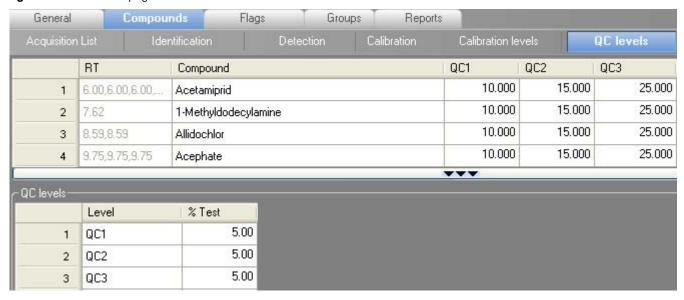


Table 35. QC levels page parameters

| Parameter | Description |
|-----------|---|
| RT | Retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| Compound | The compound name. |
| QC1-QCn | User-defined quality control levels for the compound. |
| QC levels | |
| Level | User-defined quality control level names. |
| % Test | A value for the acceptable difference (as a percentage) between the known amount and calculated (measured) amount of each QC level. |

Real Time Viewer

Use the Real Time Viewer page to specify which traces display in the real-time status pane when you perform acquisition in the Acquisition mode or when you acquire a development batch in the Method Development mode. See "Real-time Display" on page 246.

Figure 41. Real Time Viewer page

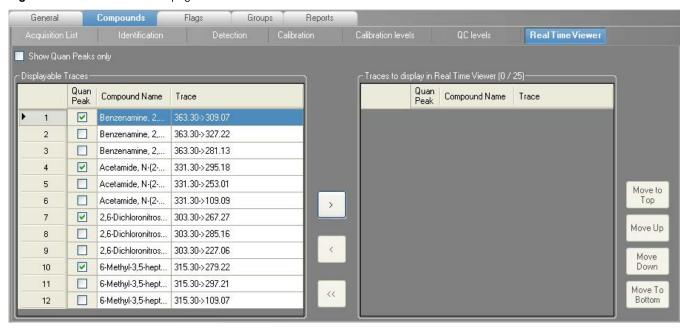


Table 36. Real Time Viewer page parameters (Sheet 1 of 2)

| Parameter | Description | |
|-------------------------|--|--|
| Show Quan Peaks Only | Displays only quan peaks in the compounds list. Quan peaks are indicated with a check in the Quan Peak column. | |
| | ✓ Show Quan Peaks only | |
| | Displayable Traces— | |
| | Quan Peak Compound Name Trace | |
| | ▶ 1 | |
| | 2 | |
| Displayable Traces | | |
| Quan Peak | Checks indicate quan peak traces. Unchecked traces indicate confirming ions. | |
| Compound Name | Names of all compounds in the method. | |
| Trace | Lists the simple mass or precursor mass for all traces—both quan peak and confirming ion—for each compound. | |
| > | Moves the selected trace to the Traces to Display in Real Time Viewer pane. | |
| < | Moves the selected trace to the Displayable Traces pane. | |

Table 36. Real Time Viewer page parameters (Sheet 2 of 2)

| Parameter | Description | |
|---|--|--|
| << | Moves all traces to the Displayable Traces pane. To move multiple traces to the Traces to Display pane, hold down the SHIFT key, select multiple traces, and then click | |
| Traces to Display in Real Time Viewer (0/25) | List the traces to be displayed and the display order in the real-time viewer in the Acquisition mode. Maximum number of traces is 25. | |
| Move to Top | Moves the selected trace to the top of the Traces to Display list and the second position in the real-time display. The TIC is always the first position in the real-time display. | |
| Move Up | Moves the selected trace up one position in the list. | |
| Move Down | Moves the selected trace down one position in the list. | |
| Move to Bottom | Moves the selected trace to the bottom of the list. | |

Using the Shortcut Menu Commands

Each page on the Compounds page (except the Acquisition List page) uses right-click shortcut menu commands to display or hide the retention column, remove compounds from the method, copy and paste data, or save the compound list to a .csv file.

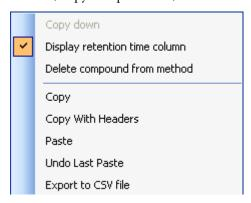


Table 37. Compounds page shortcut menu (Sheet 1 of 2)

| Command | Description |
|-------------------------------|--|
| 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. See Appendix B, "Using Copy Down and Fill Down." |
| Display retention time column | Displays or hides the RT column in the compound list. |
| Delete compound from method | Removes the selected compound from the current master method. |

4 Using the Method Development Mode

Working with Master Methods

Table 37. Compounds page shortcut menu (Sheet 2 of 2)

| Command | Description | | |
|--------------------|--|-------------------------|-----------------------------|
| Сору | Copies the data in the selected rows or columns to the Clipboard. Use this command to copy compound information to another application, such as an Excel spreadsheet. You cannot paste this data back into the method development compound list, but you can paste this copied data into other areas of the TraceFinder application. | | |
| Copy With Headers | associated column headers to the Clipboard. Use this common to copy sample information to another application, such a Excel spreadsheet. For example: Sample type | | oard. Use this command |
| | Matrix Blank | | |
| | Cal Std | | Sample type |
| | Chk Std | | Unknown/TIC |
| | Unknown/TIC Unknown/TIC | | Unknown/TIC |
| | Copy With Headers from TraceFinder | Pi | aste into Excel spreadsheet |
| | You cannot paste to compound list. | this data back into the | e method development |
| Paste | Pastes a single column of copied data from another application, such as an Excel spreadsheet, into the selected column. The pasted data must be valid data for the selected column. | | |
| Undo Last Paste | Removes the last pasted item in the method development compound list. | | |
| Export to CSV file | Opens the Save As dialog box where you can save the current compound list to a .csv file. | | |

Editing the Flags Page

Use the Flags page to set limits and ranges so the TraceFinder application can review the data and results as an aid to final approval.

From the Flags page of the Master Method View, you can access these additional pages:

- Limits
- Calibration
- Chk Std
- Matrix Blank
- Solvent Blank
- ISTD
- Surrogate
- Lab Control (Environmental and Food Safety)
- Method Val (Environmental and Food Safety)
- Matrix Spike (Environmental and Food Safety)

Limits

Use the Limits page to define levels of review for quantified results. Quantified results appear on printed and electronic reports. You can also define when a quantified value is reported instead of reporting less than a particular limit.

Figure 42. Limits page



Table 38. Limits page parameters

| Parameter | Description |
|--------------------------|--|
| RT | Retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| Compound | The compound name. |
| LOD (Detection limit) | Limit of detection. The lowest amount that can be detected. Usually derived from a method detection limit (mdl) study. |
| LOQ (Quantitation limit) | Limit of quantitation. The lowest amount that can be confidently and accurately quantitated. This is usually the lowest calibration amount. |
| LOR (Reporting limit) | Limit of reporting. Also called cutoff in some industries. This is the lowest amount that can be reported, as determined by each laboratory's standard operating practices. |
| ULOL (Linearity limit) | Upper limit of linearity. This is usually the highest calibrator amount. |
| Carryover limit | The highest amount of a substance that does not leave a residual amount in the instrument. If a substance has a carryover limit of 5, amounts higher than 5 usually dirty the instrument and leave residue behind, tainting the following sample. A carryover limit of less than 5 does not leave any residual amounts of the substance. |

Calibration

Use the Calibration page to define acceptable criteria for initial calibration. The TraceFinder application makes the evaluation by comparing the initial calibration results for each compound found in the sample to the values defined on this page.

On the Calibration report, the application flags the calculated values for internal standard compounds that exceed these limits.

Figure 43. Calibration page



Table 39. Calibration page parameters

| Parameter | Description |
|------------------|--|
| RT | Retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| Compound | The compound name. |
| R^2 threshold | The minimum correlation coefficient (r ²) for an acceptable calibration (when in linear or quadratic mode). |
| Max RSD (%) | The maximum relative standard deviation (RSD) for an acceptable calibration (when in average RF mode). |
| Min RF | The minimum average response factor (RF) for an acceptable calibration (when in average RF mode). |
| Max Amt Diff (%) | The maximum deviation between the calculated and theoretical concentrations of the calibration curve data points (when in linear or quadratic mode). |

Chk Std

Use the Chk Std page to review the calibration on an ongoing basis. The TraceFinder application makes the evaluation by comparing the quality check standard results for each compound in the sample to the initial calibration using values defined on this page.

On the Check Standard report, the TraceFinder application flags the calculated values for internal standard compounds that exceed these limits.

For linear and quadratic modes, the maximum difference for the calculated concentration in the Chk Std sample versus the theoretical value is set on the QC levels page of the Compounds page.

Figure 44. Chk Std page



Table 40. Chk Std page parameters

| Parameter | Description |
|-----------------|--|
| RT | Retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| Compound | The compound name. |
| Max RF Diff (%) | The maximum deviation between the response factor (RF) of the Chk Std sample and the average response factor from the calibration (when in average RF mode). |
| Min RF | The minimum response factor for the Chk Std sample (when in average RF mode). |

Matrix Blank

Use the Matrix Blank page to define acceptable levels of target compounds in blank samples. The TraceFinder application makes the evaluation by comparing the calculated concentration for each compound in the sample to the maximum concentration defined on this page. You can enter the maximum concentration as a percentage of a flag value or as a specified value.

On the Matrix Blank report, the application flags the calculated values for target compounds that exceed these limits.

To specify the maximum concentration as a percentage

- 1. From the Method column list, select one of the following methods:
 - None
 - Concentration
 - % of LOD
 - % of LOQ
 - % of LOR
- 2. In the Percentage column, type a percentage value.

❖ To specify the maximum concentration as an absolute value

- 1. From the Method column list, select **Concentration**.
- 2. In the Max Conc column, type an absolute value.

❖ To specify no maximum concentration

From the Method column list, select **None**.

Figure 45. Matrix Blank page



Table 41. Matrix Blank page parameters (Sheet 1 of 2)

| Parameter | Description | |
|-----------|--|--|
| RT | Retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. | |
| Compound | The compound name. | |

4 Using the Method Development Mode

Working with Master Methods

Table 41. Matrix Blank page parameters (Sheet 2 of 2)

| Parameter | Description |
|------------|--|
| Method | The evaluation process used for comparing the calculated concentration. You can specify no maximum, a specific concentration, or a percentage of the LOR, LOD, or LOQ. |
| Percentage | The percentage of the LOR, LOD, or LOQ if you are using the percentage approach. |
| Max Conc | The maximum concentration if you are using an absolute value. |

Solvent Blank

Use the Solvent Blank page to view or edit QC values for solvent reporting. The evaluation is made by comparing the calculated response for each compound in the sample to the maximum response defined on this page.

On the Solvent Blank report, the TraceFinder application flags the calculated values for target compounds that exceed these limits.

Figure 46. Solvent Blank page



Table 42. Solvent Blank page parameters

| Parameter | Description |
|-------------|--|
| RT | Retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| Compound | The compound name. |
| Method | The evaluation process to use as a response for the quan ion only (Quan Ion RT) or a summed response for the quan ion and any confirming ions (All Ion RT). To deactivate the solvent blank test for a specific compound, select None . |
| Upper Limit | Specifies an upper limit for each compound in the sample when you select an evaluation process. These values are not concentrations; they are raw response values. |

ISTD

Use the ISTD page to review the response and retention time of internal standards (if available). The TraceFinder application makes the evaluation by comparing the area and retention time results for each internal standard compound in the sample to a specified range.

If all of your target compounds are set to external calibration mode or you have not identified any compounds as internal standards, this page does not show any values.

Figure 47. ISTD page



Table 43. ISTD page parameters

| Parameter | Description |
|------------------------------------|---|
| RT | Retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| Compound | The compound name. |
| Min recovery (%) Max recovery (%) | The minimum and maximum percent recoveries for the internal standards to define an acceptable range. For check standards, the TraceFinder application compares the response of each internal standard in each sample to a range around the average of the responses of that compound in all of the calibration standards. For all other samples, the application calculates the comparison range around the check standard responses if a check standard is available in the batch. If no check standard is available, the application tests against the initial calibration. |
| Min RT (-min) | The minimum and maximum drift (in minutes) for the internal standards to define an acceptable range. For check standards, the TraceFinder application compares the retention time of each internal standard in each sample to a range around the average of the retention times of that compound in all of the calibration standards. For all other samples, the application calculates the comparison range around the check standard retention times if a check standard is available in the batch. If no check standard is available, the application tests against the initial calibration. |
| Max RT (+min) | |
| CV Test (%) | Coefficient of Variance test. |

Surrogate

Use the Surrogate page to view or edit values for surrogate recovery reporting. The application makes the evaluation by comparing the calculated concentration for each compound in the sample to the theoretical concentration and range defined on this page.

You can use these parameters to evaluate the performance of your method. For this evaluation, prepare, analyze, and evaluate a number of samples (typically 4 to 10) to document method accuracy and precision as a comprehensive whole.

On the Surrogate Recovery report, the application flags the calculated values for method validation compounds that exceed these limits.

Figure 48. Surrogate page



Table 44. Surrogate page parameters

| Parameter | Description |
|------------------|---|
| RT | Retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| Compound | The compound name. |
| Theo Conc | Values for the compounds that represent the expected theoretical concentration of that compound in the sample. |
| Min recovery (%) | A range of the allowable minimum recovery percentage and the maximum recovery percentage that the application can determine by comparing the observed calculated concentration in the analysis to the expected concentration. Each method validation compound can have its own values for these fields, independent of other method validation compounds. |
| Max recovery (%) | |
| Max RSD (%) | The maximum relative standard deviation of the set of observed concentrations for a component across the set of method validation samples. |

Lab Control (Environmental and Food Safety)

Use the Lab Control page to view and edit QC values for lab control sample (LCS) and lab control sample duplicate (LCSD) analyses. The application makes the evaluation by comparing the calculated concentration for each compound in the sample to the theoretical concentration and range defined on this page.

You can prepare samples (typically known as clean matrices) as LCS or LCSD. These represent samples where you have added known concentrations of target analytes. To define an LCS and its duplicate in a batch, select the appropriate sample type and a common sample ID.

On the Lab Control report, the application flags the calculated values for spiked compounds that exceed these limits.

Figure 49. Lab Control page



Table 45. Lab Control page parameters

| Parameter | Description |
|------------------------------------|---|
| RT | Retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| Compound | The compound name. |
| Theo Conc | Values for the lab control compounds that represent the expected theoretical concentration of that compound in the sample. |
| Min recovery (%) Max recovery (%) | A range of the allowable minimum recovery percentage and the maximum recovery percentage that the application can determine by comparing the observed calculated concentration in the analysis to the expected concentration. Each LCS or LCSD compound can have its own values for these fields, independent of other LCS or LCSD compounds. |
| Max RPD | Specifies a maximum value for relative percent difference (RPD) between two spiked samples. |

Method Val (Environmental and Food Safety)

Use the Method Val page to view or edit QC values for method validation reporting. The application makes the evaluation by comparing the calculated concentration for each compound in the sample to the theoretical concentration and range defined on this page.

You can use these parameters to evaluate the performance of your method. For this evaluation, prepare, analyze, and evaluate a number of samples (typically 4 to 10) to document method accuracy and precision as a comprehensive whole. To define a method validation sample in the batch, select the appropriate sample type.

On the Method Validation report, the application flags the calculated values for method validation compounds that exceed these limits.

Figure 50. Method Val page



Table 46. Method Val page parameters

| Parameter | Description |
|------------------|--|
| RT | Retention time for the compound. |
| Compound | The compound name. |
| Theo Conc | Values for the compounds that represent the expected theoretical concentration of that compound in the sample. |
| Min recovery (%) | A range of the allowable minimum recovery percentage and the |
| Max recovery (%) | maximum recovery percentage that the application can determine by comparing the observed calculated concentration in the analysis to the expected concentration. Each method validation compound can have its own values for these fields, independent of other method validation compounds. |
| Max RSD (%) | The maximum relative standard deviation of the set of observed concentrations for a component across the set of method validation samples. |

Matrix Spike (Environmental and Food Safety)

Use the Matrix Spike page to view or edit QC values for matrix spike and matrix spike duplicate analyses. The application makes the evaluation by comparing the calculated concentration for each compound in the sample (after subtracting the original sample value) to the theoretical concentration and range defined on this page.

To evaluate matrix spike or matrix spike duplicate compounds, prepare samples as MS or MSD. These represent samples where you have added known concentrations of target analytes. To define a sample, its MS, and its MSD in the batch, select the appropriate Sample Type and a Sample ID.

Sample IDs must be unique. Duplicating Sample IDs can cause incorrect samples to be included in reports.

On the MS/MSD report, the application flags the calculated values for spiked compounds that exceed these limits.

Figure 51. Matrix Spike page



Table 47. Matrix Spike page parameters

| Parameter | Description |
|------------------|--|
| RT | Retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| Compound | The compound name. |
| Theo Conc | Values for the matrix spike compounds that represent the expected theoretical concentration of that compound in the sample. You can apply or not apply the dilution factor for a sample to the calculated matrix spike concentrations. |
| Min recovery (%) | A range of the allowable minimum recovery percentage and the |
| Max recovery (%) | maximum recovery percentage that can be determined by comparing the observed calculated concentration in the analysis to the expected concentration. Each matrix spike sample can have its own values for these fields, independent of other matrix spike samples. |
| Max RPD | Specify a maximum value for relative percent difference (RPD) between two spiked samples. |

Editing the Groups Page

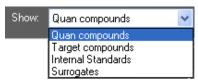
Use the Groups page of the Master Method View to organize compounds into functional or logical groups. You can use these groups for creating a subset of target compounds. For a detailed description of all the features on the Groups page, see "Groups page parameters" on page 172.

For quantitative processing, the TraceFinder application processes all compounds in the method and stores the complete result set, but only those in the selected group are visible in the Acquisition mode. Limiting the displayed compounds to those in the selected group can be useful when working with a master method containing a large list of compounds, only some of which are required for analysis in certain samples. In that case, the application requires only a single method and can reduce the results. To display only those compounds to be used in quantitative processing, select **Quan Compounds** from the Show list.

You can create multiple groups and include the same compound in more than one group.

❖ To create a group

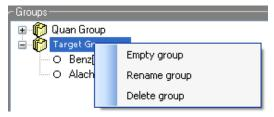
1. From the Show list, select the type of compounds you want to view.



2. At the bottom of the Groups area, click **Add Group**.

The Add a New Group dialog box opens.

- Type a name for the new group and click **OK**.The new group appears in the Groups area.
- 4. Drag a compound from the Compounds area onto a group name (as if you were moving files into a folder).
- 5. To remove all the compounds from a group, rename the group, or delete it, right-click the group name and choose from the shortcut menu.



6. To remove a single compound, right-click the compound name in the group and choose **Remove from Group** from the shortcut menu.



Working with Master Methods

Figure 52. Groups page

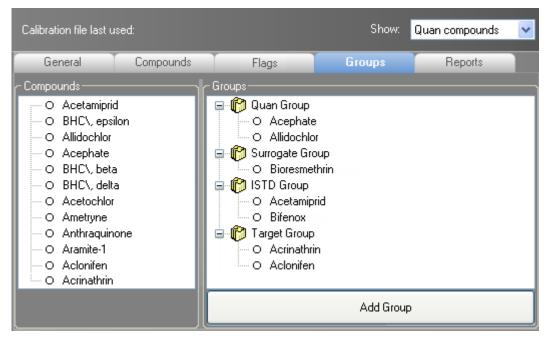


Table 48. Groups page parameters

| Parameter | Description |
|-------------------|--|
| Compounds | Lists all available compounds. |
| Groups | Lists all available groups. |
| Add Group | Opens the Add a New Group dialog box where you can create a new group. |
| Shortcut menu | |
| Empty group | Removes all compounds from the selected group. |
| Rename group | Changes the name of the selected group. |
| Delete group | Removes the selected group and all the compounds in it. |
| Remove from group | Removes the selected compound from its group. |

Editing the Reports Page

Use the Reports page to specify how you want to save or print your reports. For a detailed description of the features on the Reports page, see "Reports page parameters" on page 176.

For the Quantitation Report type, you can modify quan report, user interface, quan flag, and surrogate correction options in the Report Options pane.

This section includes instructions for the following tasks:

- Specifying Report Formats
- Specifying Quan Report Settings
- Specifying Target Screening Settings

Specifying Report Formats

- For each standard report, you can create a hardcopy printout, a PDF file, or an XML file.
- For each custom report, you can create a hardcopy printout or an Excel Macro-Enabled Workbook (.xlsm) file.
- For each target screening report, you can create a hardcopy printout or a PDF file.

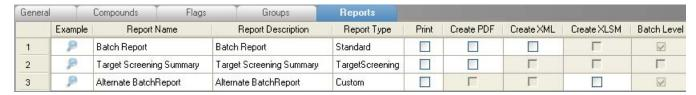
In addition to the report type, you can specify a report description for each of your reports. The default report description is the report name.

To specify standard report types and output formats

1. Click the **Reports** tab.

The Reports page displays the following columns for all configured reports:

- The Example icon that opens an example PDF of the report type
- Report Name, Report Description, and Report Type
- Options to create a hardcopy, PDF file, or XML file for standard report types
- Options to create a hardcopy or Excel Macro-Enabled Workbook file for custom report types
- Options to create a hardcopy or PDF file for target screening report types
- The Batch Level option that indicates which reports are batch-level reports



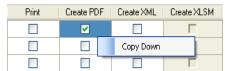
For information about configuring which reports are available when you create a master method or which reports create a batch-level report, see "Specifying the Reports Configuration" on page 58.

Working with Master Methods

2. To edit the Report Description, double-click the name and type your new name.

The TraceFinder application uses this name for all reports that use this master method. You cannot edit the Report Description from other report views.

- 3. To specify the type of report output to create for each report type, select the check box in the appropriate column.
- 4. To duplicate the output type for all reports, click the cell to select it, then right-click and choose **Copy Down** from the shortcut menu.



All check boxes in the column below the selected cell duplicate the selected or cleared state of the selected cell. This action applies only to reports where this output format is available.

By default, all report types are cleared.

Figure 53. Reports page

| | Example | Report Name | Report Description | Report Type | Print | Create PDF | Create XML | Create XLSM | Batch Level |
|----|---------|--|-----------------------------|-----------------|-------|------------|------------|-------------|-------------|
| 1 | P | Batch Report | Batch Report | Standard | | | | | ✓ |
| 2 | P | Batch Report Rev 1 | Batch Report Rev 1 | Standard | | | | | ✓ |
| 3 | P | Blank Report | Blank Report | Standard | | | | | |
| 4 | P | Calibration Report | Calibration Report | Standard | | | | | ✓ |
| 5 | P | Check Standard Report | Check Standard Report | Standard | | | | | Г |
| 6 | P | Chromatogram Report | Chromatogram Report | Standard | | | | | Г |
| 7 | P | Compound Calibration Report | Compound Calibration Re | Standard | | | | | Г |
| 8 | P | Compound Calibration Report - Alternate | Compound Calibration Re | Standard | | | | | Г |
| 9 | P | Confirmation Report | Confirmation Report | Standard | | | | Г | Г |
| 10 | P | Confirmation Report 2 | Confirmation Report 2 | Standard | | | | | Г |
| 11 | P | High Density Calibration Report | High Density Calibration R | Standard | | | | Г | V |
| 12 | P | High Density Internal Standard Report | High Density Internal Stan | Standard | | | | | Г |
| 13 | P | High Density Internal Standard Report Long | High Density Internal Stan | Standard | | | | | Г |
| 14 | P | High Density Sample Report 1 | High Density Sample Rep | Standard | | | | | Г |
| 15 | P | High Density Sample Report 1 Long | High Density Sample Rep | Standard | | | | Г | Г |
| 16 | P | High Density Sample Report 2 | High Density Sample Rep | Standard | | | | Г | Г |
| 17 | P | High Density Sample Report 2 Long | High Density Sample Rep | Standard | | | | | Г |
| 18 | P | High Density Sample Report 3 | High Density Sample Rep | Standard | | | | Г | Г |
| 19 | P | High Density Sample Report 3 Long | High Density Sample Rep | Standard | | | | | Г |
| 20 | P | Internal Standard Summary Report | Internal Standard Summar | Standard | | | | Г | Г |
| 21 | P | Ion Ratio Failure Report | Ion Ratio Failure Report | Standard | | | | | Г |
| 22 | P | Manual Integration Report | Manual Integration Report | Standard | | | | | Г |
| 23 | P | Method Report | Method Report | Standard | | | | | V |
| 24 | P | Quantitation Report | Quantitation Report | Standard | | | | | Г |
| 25 | P | Quantitation Report - 2 | Quantitation Report - 2 | Standard | | | | | Г |
| 26 | P | Solvent Blank Report | Solvent Blank Report | Standard | | | | Г | Г |
| 27 | P | Surrogate Recovery Report | Surrogate Recovery Report | Standard | | | | | Г |
| 28 | P | TIC Report | TIC Report | Standard | | | | Г | Г |
| 29 | P | TIC Summary Report | TIC Summary Report | Standard | | | | | Г |
| 30 | P | Target Screening Long Report | Target Screening Long R | TargetScreening | | | Г | Г | Г |
| 31 | P | Target Screening Summary Report | Target Screening Summar | TargetScreening | | | Г | Г | Г |
| 32 | P | AltCalibrationReport | AltCalibrationReport | Custom | | Г | Г | | V |
| 33 | P | Alternate BatchReport | Alternate BatchReport | Custom | | Г | Г | | V |
| 34 | P | Alternate CalibrationReport | Alternate CalibrationReport | Custom | | Г | | | ✓ |
| 35 | P | Alternate ConfirmationReport | Alternate ConfirmationRep | Custom | | Г | | | Г |
| 36 | P | Alternate MatrixSpikeReport | Alternate MatrixSpikeReport | Custom | | | Г | | Г |

4 Using the Method Development Mode Working with Master Methods

Figure 54. Reports page, continued

| P | Alternate SampleReport | Alternate SampleReport | Custom | | Г | Г | | Г |
|---|------------------------------|---|---|--|--|--|-------------------------|--|
| P | Alternate SummaryReport | Alternate SummaryReport | Custom | | Г | Г | | V |
| P | BatchReport | BatchReport | Custom | | Г | Г | | V |
| P | BlankReport | BlankReport | Custom | | Г | Г | | Г |
| P | CalibrationDensityReport | CalibrationDensityReport | Custom | | Г | Г | | ~ |
| P | CalibrationReport | CalibrationReport | Custom | | Г | Г | | ~ |
| P | CheckStandardReport | CheckStandardReport | Custom | | Г | Г | | Г |
| P | CompoundCalibrationReport | CompoundCalibrationReport | Custom | | | | | |
| P | ConfirmationReport | ConfirmationReport | Custom | | | | | |
| P | ConfirmationReport2 | ConfirmationReport2 | Custom | | | | | |
| P | HighDensitySampleReport1Long | HighDensitySampleRepor | Custom | | | | | |
| P | HighDensitySampleReport2Long | HighDensitySampleRepor | Custom | | | | | |
| P | HighDensitySampleReport3Long | HighDensitySampleRepor | Custom | | | | | |
| P | HighDensitySampleReport4 | HighDensitySampleReport4 | Custom | | | | | |
| P | HighDensitySampleReport5 | HighDensitySampleReport5 | Custom | | Г | Г | | Г |
| P | QuantitationReport | QuantitationReport | Custom | | Г | Г | | |
| P | SteroidAnalysisReport | SteroidAnalysisReport | Custom | | Г | Г | | Г |
| | | Alternate SummaryReport BatchReport BlankReport CalibrationDensityReport CalibrationReport CheckStandardReport CompoundCalibrationReport ConfirmationReport ConfirmationReport2 HighDensitySampleReport2Long HighDensitySampleReport3Long HighDensitySampleReport4 HighDensitySampleReport4 HighDensitySampleReport5 QuantitationReport | Alternate SummaryReport BatchReport BlankReport BlankReport CalibrationDensityReport CalibrationDensityReport CalibrationReport CalibrationReport CalibrationReport CheckStandardReport CompoundCalibrationReport ConfirmationReport ConfirmationReport ConfirmationReport ConfirmationReport ConfirmationReport ConfirmationReport ConfirmationReport HighDensitySampleReport1Long HighDensitySampleReport3Long HighDensitySampleReport3Long HighDensitySampleReport4 HighDensitySampleReport4 HighDensitySampleReport4 HighDensitySampleReport5 QuantitationReport QuantitationReport QuantitationReport | Alternate SummaryReport Alternate SummaryReport BatchReport BlankReport Custom CalibrationDensityReport CalibrationReport CalibrationReport CalibrationReport CalibrationReport CalibrationReport CalibrationReport CheckStandardReport CompoundCalibrationReport ConfirmationReport ConfirmationReport ConfirmationReport ConfirmationReport ConfirmationReport ConfirmationReport ConfirmationReport Custom ConfirmationReport Custom HighDensitySampleReport1Long HighDensitySampleReport. Custom HighDensitySampleReport3Long HighDensitySampleReport. HighDensitySampleReport4 HighDensitySampleReport4 HighDensitySampleReport5 Custom Custom HighDensitySampleReport4 Custom HighDensitySampleReport4 Custom Custom Custom HighDensitySampleReport5 HighDensitySampleReport5 Custom Custom Custom Custom Custom Custom HighDensitySampleReport4 Custom Custom | Alternate SummaryReport Alternate SummaryReport BatchReport BatchReport BlankReport Custom CalibrationDensityReport CalibrationPensityReport CalibrationPensityReport CalibrationPenort CalibrationPenort CalibrationPenort CalibrationPenort CalibrationPenort CalibrationPenort Custom CheckStandardReport CompoundCalibrationReport CompoundCalibrationReport ConfirmationReport ConfirmationReport ConfirmationReport ConfirmationReport ConfirmationReport ConfirmationReport Custom HighDensitySampleReport1Long HighDensitySampleReport Custom HighDensitySampleReport2Long HighDensitySampleReport Custom HighDensitySampleReport Custom HighDensitySampleReport Custom HighDensitySampleReport Custom HighDensitySampleReport Custom HighDensitySampleReport Custom Custom HighDensitySampleReport Custom Custom HighDensitySampleReport Custom Custom HighDensitySampleReport Custom QuantitationReport Custom QuantitationReport Custom | Alternate SummaryReport Alternate SummaryReport BatchReport BatchReport BlankReport Custom CalibrationDensityReport CalibrationDensityReport CalibrationBeport CalibrationReport CalibrationReport CalibrationReport CalibrationReport CheckStandardReport CompoundCalibrationReport ConfirmationReport Custom ConfirmationReport Custom ConfirmationReport Custom ConfirmationReport Custom ConfirmationReport Custom Cu | Alternate SummaryReport | Alternate SummaryReport Alternate SummaryReport Custom BatchReport BatchReport BlankReport BlankReport Custom CalibrationDensityReport Custom CalibrationReport CalibrationReport Custom CheckStandardReport Custom CompoundCalibrationReport CompoundCalibrationReport ConfirmationReport ConfirmationReport ConfirmationReport ConfirmationReport ConfirmationReport ConfirmationReport Custom |

 Table 49.
 Reports page parameters

| Parameter | Description |
|--------------------|---|
| Example | Opens a PDF that displays an example of the report type. |
| Report Name | The name of a report. |
| Report Description | The user-defined description to be used on a report. |
| Print | Sends reports to the printer. |
| Create PDF | Saves reports as PDF files. Available only for standard reports. |
| Create XML | Exports reports in XML format. Available only for standard reports. |
| Create XLSM | Exports reports in Excel Macro-Enabled Workbook (.xlsm) format. Available only for custom reports. |
| Batch Level | Rather than creating separate reports for each sample, the application uses a composite of the data from all the samples to create a single report for the entire batch. Batch-level reports are prepended with a B to differentiate them. You cannot select this option from the Reports page. You must select the Batch Level option for the report in the report configuration. See "Specifying the Reports Configuration" on page 58. |

Specifying Quan Report Settings

Use the options on the Quan Report Settings page to choose parameters for flagging values and displaying information in standard report types.

Use the following procedures:

- To specify quantitation limits
- To specify user interface options
- To specify quantitation flag options
- To correct surrogates
- To track the use of the tune file

❖ To specify quantitation limits



 To report the calculated concentration at all times or only when the quantified value exceeds LOD, LOQ, or LOR, choose the appropriate value from the Report Concentration list.

For a description of concentration limits, see "Limits" on page 160.

- 2. To select the number of decimal places to report for calculated concentrations, set the value in the Decimal Places to be Reported box.
- 3. To include a chromatogram of the sample in the Quantitation Report, select the **Show Chromatogram on Quantitation Report** check box.
- 4. To display only valid compounds, select the **Display Compounds Above Set Limit** check box.

❖ To specify user interface options



- 1. To shade a compound row on any of the reports if a value fails one of the criteria used for evaluation, select the **Shade Row when Sample is Outside of Evaluation Criteria** check box.
- 2. To separate the ion overlay pane from the confirming ion plots, select the **Separate Ion Overlay Display** check box.
- 3. To use an alternate format for the Calibration Report designed to print more concisely and limit the report to a maximum of 7 calibration standards, select the **Use Alternate Calibration Report Format** check box.
- 4. To display flags and a legend on high density reports, select the **Display Quan Flags and Legend** check box.

To specify quantitation flag options



Select the values you want to display in the report.

Values are above or below the limits defined on the Quan page.

These flags appear on a variety of reports and are defined in the "Quan Report Settings page parameters" on page 180.

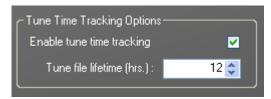
To correct surrogates



Select the Correct Surrogates check box.

The TraceFinder application applies the conversion factor (specified in the sample row in the batch) to the sample's calculated concentrations for surrogates as the conversion factor is applied to target compounds.

❖ To track the use of the tune file



1. Select the **Enable Tune Time Tracking** check box.

This option tracks the number of hours between the last instrument tune and each sample acquisition.

2. In the Tune File Lifetime box, enter the number of hours you want to allow between the last instrument tune and a sample acquisition.

Any sample acquired outside this maximum allowable time is flagged on the Batch report.

Working with Master Methods

Figure 55. Quan Report Settings page

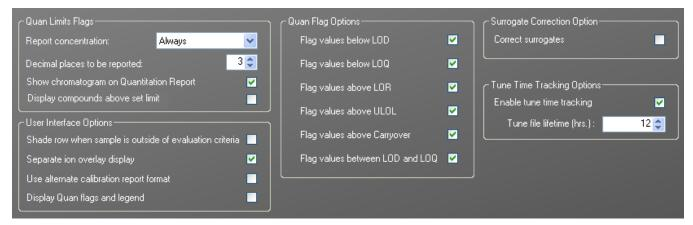


Table 50. Quan Report Settings page parameters (Sheet 1 of 2)

| Parameter | Description |
|---|--|
| Quan Limits Flags | |
| Report concentration | Reports the concentration at all times or only when the quantified value exceeds either the limit of detection (LOD), the limit of quantitation (LOQ), or the limit of reporting (LOR). Report concentration: Always, >LOD, >LOQ, or >LOR. |
| Decimal places to be reported | Number of decimal places to be included in the report. Maximum value is 6. |
| Show chromatogram on Quantitation Report | Displays a chromatogram (TIC trace) of the sample on the quantitation report. |
| Display compounds above set limit | Prints only the positive compounds in a sample. If a compound is above the specified Quan Flag Options limits, the TraceFinder application reports the compound. |
| User Interface Options | |
| Shade row when sample is outside of evaluation criteria | Shades a compound row on any of the reports if a value fails one of the criteria used for evaluation. |
| Separate ion overlay display | Separates the ion overlay pane from the confirming ion plots in data review. |
| Use alternate calibration report format | Uses an alternate format for the Calibration Report that is designed to print more concisely (this report is limited to a maximum of 7 calibration standards). |
| Display Quan flags and legend | Displays manual flags, confirming manual flags, quan flags, and a legend on high density reports. |

Table 50. Quan Report Settings page parameters (Sheet 2 of 2)

| Parameter | Description | | | | | |
|------------------------------------|---|--|--|--|--|--|
| Quan Flag Options | Values that are above or below limits defined on the Limits page. These flags appear on a variety of reports. | | | | | |
| Flag values below LOD | Flags values below the limit of detection (LOD). | | | | | |
| Flag values below LOQ | Flags values below the limit of quantitation (LOQ). | | | | | |
| Flag values above LOR | Flags values above the limit of reporting (LOR). | | | | | |
| Flag values above ULOL | Flags values above the upper limit of linearity (ULOL). | | | | | |
| Flag values above Carryover | Flags values above the carryover limit. | | | | | |
| Flag values between LOD and LOQ | Flags values between the limit of detection and the limit of quantitation known as the J flag. | | | | | |
| Surrogate Correction Opti | ion | | | | | |
| Correct surrogates | Applies the conversion factor (specified in the sample row in the batch) to the sample's calculated concentrations for surrogates as the conversion factor is applied to target compounds. For example, if you added surrogates to the sample as part of sample preparation and you require a dilution for analysis, the TraceFinder application dilutes the surrogates and target compounds and applies a dilution correction to correct for this dilution. However, if you added surrogates after a dilution has occurred, then you can leave the option cleared so that, while the target compounds are corrected for the dilution, the surrogates are reported "as is." | | | | | |
| Tune Time Tracking Optio | Tune Time Tracking Options | | | | | |
| Enable tune time tracking | Tracks the number of hours between the last instrument tune and each sample acquisition. | | | | | |
| Tune file lifetime | Specifies the maximum number of hours between the last instrument tune and a sample acquisition. Any sample acquired outside this maximum allowable time is flagged on the Batch report. | | | | | |

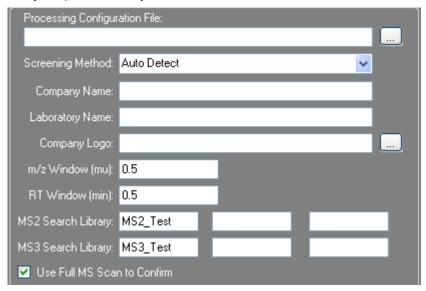
Specifying Target Screening Settings

Use the options on the Target Screening Settings page to set the parameters required to produce Target Screening reports. For a detailed description of the features on the Target Screening Settings page, see "Target Screening Settings page parameters" on page 185.

The TraceFinder application use these parameters to process a raw data file and create a report similar to a ToxID report. See "Example Target Screening Summary Report" on page 187.

Use the following procedures:

- To specify the default parameters
- To calculate and report semi-quantitative results
- To specify the ion ratio calculation method
- To specify the exact mass window
- To specify the Exactive parameters
- ❖ To specify the default parameters



- 1. Click the Processing Configuration File browse button and select a configuration file (.csv).
- 2. From the Screening Method list, select one of these compound screening methods.
 - •(Default) Auto Detect
 - •Based on Full MS2 scans
 - •Based on SRM and MS2 scans
 - •Based on MS2 and MS3 scans
 - •Based on MS3 scans
 - •Based on accurate mass scans
 - •Based on SRM scans
 - •Based on Exactive screening method

- 3. Type the name of the company to print on the report.
- 4. Type the name of the laboratory to print on the report.
- 5. Click the Company Logo browse button and select a graphic file (.jpg, .gif, or .bmp) to print on the report.
- 6. In the *m*/*z* Window box, enter a value for the window above and below the m/z value for the compounds.
- 7. In the RT Window box, enter a value for the window above and below the retention time value for the compounds.
- 8. In the MS2 Search Library boxes, type the names of as many as three search libraries for searching MS/MS spectra.
- 9. In the MS3 Search Library boxes, type the names of as many as three search libraries for searching MS³ spectra.
- 10. Select the **Use Full MS Scan to Confirm** check box if you want to confirm library search results with parent ion peak detection in the full scan.

When the application does not detect a peak in the full scan, the compound is not reported as a hit.

❖ To calculate and report semi-quantitative results



- 1. In the Semi Quantitative area, do the following:
 - a. Select the **Report Semi-Quantitative Result** check box.
 - b. Type the measurement units.

The measurement units are used only for labeling purposes.

- 2. Select either the **Scan Intensity** or **Peak Area** option.
 - Scan Intensity: The application measures the intensity of the MS/MS peak without performing background subtraction.
 - Peak Area: The application measures the peak area of the reconstructed full scan chromatogram peak of the parent ion. When you select Peak Area, the Use Full MS Scan to Confirm check box is automatically selected.

To specify the ion ratio calculation method



- 1. Select the Use Scan at Peak Apex or Use Average Scan option.
 - Use Scan at Peak Apex: The application calculates the ion ratio based on the peak apex scan spectrum.
 - Use Average Scan: The application calculates the ion ratio based on the average scan spectrum over the range of the peak's half height. The relative intensity in the configuration file must use the selected scan method.
- 2. In the Ion Ratio Window (%) box, type the acceptable percentage of the intensity of the qualifier ion to the quan ion.

For example, when the Ion Ratio Window is 20% and the quan ion has an intensity/height of 100, the specified confirming ion/mass must have a height of at least 80 to be considered found.

❖ To specify the exact mass window



Type a total window width value in parts per million for the Exact Mass Window.

For example, when you expect a mass of 50 with a window of 2, the algorithm creates an XIC based on the responses of all masses from 49 to 51.

❖ To specify the Exactive parameters



1. Type values for Adduct 1, Adduct 2, and Adduct 3.

These values identify the adducts listed in and applied throughout the configuration file. Adducts are polarity sensitive. For negative ionization, enter negative adduct values.

These values default to H+, NH4+, and Na+, respectively.

- 2. To search the entire raw data file for the specified peak, do the following:
 - a. Select the **No Specified Retention Time** check box.
 - b. Select either the First Peak or Highest Peak option.

When the search finds more than one m/z match in the raw data file, the application uses the specified peak for processing.

3. Select the **Report All Compounds Listed in Configuration File** check box to report all compounds in the configuration file whether or not matches are found for them.

The default reports on only those compounds where matches are found in the raw data file. This option applies to the Exactive experiment only.

Figure 56. Target Screening Settings page

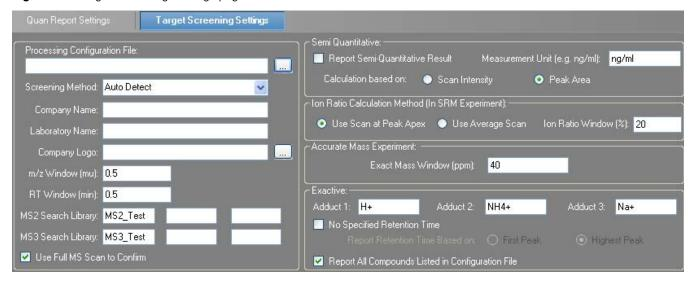


Table 51. Target Screening Settings page parameters (Sheet 1 of 2)

| Parameter | Description |
|----------------------------------|--|
| Processing Configuration File | Specifies a configuration file (.csv). |
| Screening Method | Specifies one of the following screening methods: • (Default) Auto Detect • Based on Full MS2 scans • Based on SRM and MS2 scans • Based on MS2 and MS3 scans • Based on MS3 scans • Based on accurate mass scans • Based on SRM scans • Based on Exactive screening method Note Using the Auto Detect method, the ToxID application can identify the screening experiment implemented in the acquired data file. |
| Company Name | Specifies the name of the company to print on the report. |
| Laboratory Name | Specifies the name of the laboratory to print on the report. |
| Company Logo | Specifies a graphic file (.jpg, .gif, or .bmp) to print on the report. |
| m/z Window (mu) | Specifies a value for the window above and below the m/z value for the compounds. |
| RT window (min) | Specifies a value for the window above and below the retention time value for the compounds. |

4 Using the Method Development Mode Working with Master Methods

Table 51. Target Screening Settings page parameters (Sheet 2 of 2)

| Parameter | Description |
|---|---|
| MS2 Search Library | Specifies the names of as many as three search libraries for searching MS/MS spectra. |
| MS3 Search Library | Specifies the names of as many as three search libraries for searching MS ³ spectra. |
| Use Full MS Scan to Confirm | Specifies that the application confirms library search results with parent ion peak detection in the full scan. When the application does not detect a peak in the full scan, the compound is not reported as a hit. |
| Semi Quantitative | |
| Report Semi-Quantitative Result | |
| Measurement Unit | |
| Calculation based on | Specifies one of the following calculation methods: Scan Intensity: Specifies that the application measures the intensity of the MS/MS peak without performing background subtraction. Peak Area: Specifies that the application measures the peak area of the reconstructed full scan chromatogram peak of the parent ion. When you select Peak Area, the Use Full MS Scan to Confirm check box is automatically selected. |
| Ion Ratio Calculation Me | thod (In SRM Experiment) |
| Use Scan at Peak Apex | Specifies that the application calculates the ion ratio based on the peak apex scan spectrum. |
| Use Average Scan | Specifies that the application calculates the ion ratio based on the average scan spectrum over the range of the peak's half height. The relative intensity in the configuration file must use the selected scan method. |
| Ion Ratio Window(%) | |
| Accurate Mass Experime | ent |
| Exact Mass Window | Specifies a value in parts per million for the accurate mass experiment. |
| Exactive | |
| Adduct 1-n | Specifies the adducts listed in and applied throughout the configuration file. Adducts are polarity sensitive. For negative ionization, enter negative adduct values. Defaults: Adduct 1: H+, Adduct 2: NH4+, and Adduct 3: Na+ |
| No Specified Retention Time | Specifies either First Peak or Highest Peak to use for processing when the search finds more than one m/z match in the raw data file. |
| Report All Compounds Listed in Configuration File | Specifies that in an Exactive experiment, the application reports all compounds in the configuration file whether or not matches are found for them. Default: Reports only those compounds where matches are found in the raw data file. |

Figure 57. Example Target Screening Summary Report

Your Company Name Summary Report

Raw File Name: C:\Xcalibur\examples\ToxID\Exact_Mass\Exact_Mass_Test.RAW

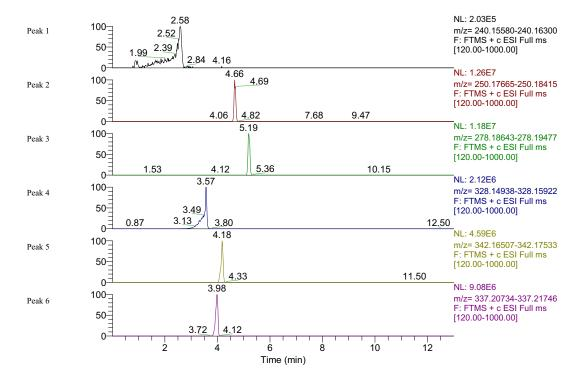
Config File Name: C:\Xcalibur\examples\ToxID\Exact_Mass\ConfigFile_Exact_mass.csv

Sample Name:

Laboratory: Your Lab Name

Acquistion Start Time: 3/24/2008 4:46:43 PM

Screening Conditions: Based on accurate mass scans. Exact mass window (ppm): 30, RT window(min): 0.50.



| Peak Number | Compound Name | Expected m/z | Detected m/z | Delta (mDa) | Delta (ppm) | Expected A | Actual RT | Intensity |
|----------------|------------------|-----------------|-----------------|----------------|----------------|------------|-----------|-----------|
| 1 | Albuterol | 240.15940 | 240.15939 | -0.0 | -0.0 | 2.58 | 2.58 | 199505 |
| 2 | Alprenolol | 250.18040 | 250.18039 | -0.0 | -0.0 | 4.50 | 4.66 | 12604499 |
| 3 | Amitriptyline | 278.19060 | 278.19061 | 0.0 | 0.0 | 5.00 | 5.19 | 11769755 |
| 4 | 6-Acetylmorphine | 328.15430 | 328.15433 | 0.0 | 0.1 | 3.30 | 3.57 | 2112090 |
| 5 | 6-Acetylcodeine | 342.17020 | 342.17035 | 0.1 | 0.4 | 4.10 | 4.18 | 4593306 |
| 6 | Acebutolol | 337.21240 | 337.21246 | 0.1 | 0.2 | 3.80 | 3.98 | 9077282 |

Creating a Method Template

In the TraceFinder application, you can create a processing method using a method template that contains the basic settings.

Use the following procedures:

- To create a method template
- To specify peak criteria
- To identify the peaks
- To specify confirming ions
- To calibrate the compounds
- To enter a note for the method
- To save the method template
- To create a method template
- 1. Click **Method Development** from the dashboard or the navigation pane.



The Method Development navigation pane opens.

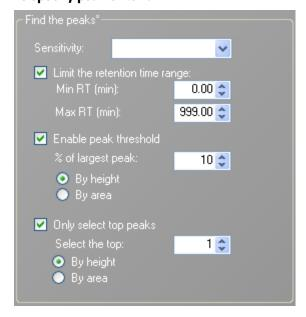
2. Click **Method View** in the navigation pane.



3. From the main menu, choose **File > New > Method Template**.

The Method Template Editor opens. For a complete description of the Method Template Editor, see "Method Template Editor dialog box" on page 195.

❖ To specify peak criteria



1. In the Find the Peaks area, select a sensitivity level.

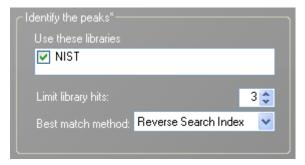
In selecting the degree of sensitivity, you define how extensively the peak detector algorithm searches for low-level peaks.

- The Genesis peak detection algorithm is provided for backward compatibility with Xcalibur 1.0 studies.
- The ICIS peak detection algorithm is designed for MS data and has superior peak detection efficiency at low MS signal levels.
- The Avalon peak detection algorithm is designed for integrating UV/Vis and analog chromatograms.
- 2. To look for peaks only in a certain range of the entire chromatogram, select the **Limit the Retention Time Range** check box and specify a retention time (RT) range.
- 3. To indicate whether to select peaks by relative height or area and the percentage of the highest peak that results in compound selection, select the **Enable Peak Threshold** check box.

To consider a peak for a processing method, the TraceFinder application uses the Enable Peak Threshold filter to determine which peaks meet the specified percentage of the largest peak.

4. To display a specific number of the largest peaks by height or area, select the **Only Select Top Peaks** check box and enter the number of peaks to display.

To identify the peaks

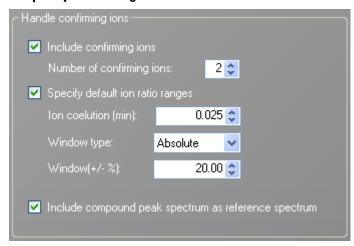


1. In the Use these Libraries box, select the libraries you want to search.

All libraries loaded on your instrument are displayed in the Use these Libraries box.

- 2. To limit the number of hits returned when the system searches a spectrum against the selected libraries, set a value in the Limit Library Hits box.
- 3. To specify how to sort the library searches, select a value from the Best Match Method list.

❖ To specify confirming ions



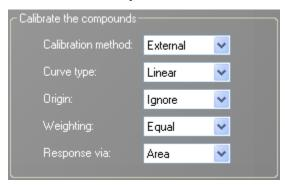
1. To set the number of confirming ions, select the **Include Confirming Ions** check box and enter a value in the Number of Confirming Ions box.

This value is the number of other ions in the spectrum whose ratio is compared to the quantitation ion. Using this ratio, you can then determine if it is the target compound or something else. This value defaults to **2** because you typically perform a 3-ion experiment with one quan mass and two confirming ions.

The system selects the most intense ion to use as the quantitation mass and uses this mass for the mathematical operations.

- 2. To define the criteria for evaluating confirming or qualifying ions, select the **Specify Default Ion Ratio Ranges** check box and set the following values:
 - a. To specify the maximum difference in retention time between a confirming ion peak and the quantification ion peak, set a value in the Ion Coelution (min) box.
 - b. To specify an absolute or relative calculation approach for determining the acceptable ion ratio range, select **Absolute** or **Relative** from the Window Type list.
 - c. To specify the acceptable ion ratio range, set a value in the Window (+/-%) box.
- 3. To include the peak spectrum in the processing method, select the **Include Compound Peak Spectrum as Reference Spectrum** check box.

To calibrate the compounds

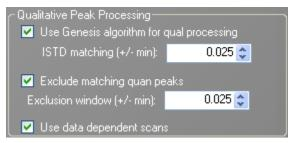


- 1. From the Calibration Method list, select Internal or External.
- 2. From the Curve Type list, select one of the following:
 - Linear: All other settings are available with this exception: When you select Include in the Origin list, all weighting values are unavailable except for Equal.
 - Quadratic: All other settings are available with this exception: When you select Include in the Origin list, all weighting values are unavailable except for Equal.
 - Average RF: No selections in the Weighting or Origin lists are available. The Weighting list is set to Equal, and the Origin list is set to Ignore.
- 3. From the Origin list, select one of the following:
 - Ignore: Specifies that the origin is not included as a valid point in the calibration curve when the curve is generated. When you select Ignore, the calibration curve might or might not pass through the origin.
 - Force: Specifies that the calibration curve passes through the origin of the data point plot when the calibration curve is generated.
 - Include: Specifies that the origin is included as a single data point in the calculation of the calibration curve. When you select Include, the calibration curve might or might not pass through the origin.

Working with Master Methods

- 4. From the Weighting list, select one of the following:
 - Equal: Specifies that the origin is included as a single data point in the calculation of the calibration curve. When you select Equal, the calibration curve might or might not pass through the origin.
 - 1/X: Specifies a weighting of 1/X for all calibration data points during the least-squares regression calculation of the calibration curve. Calibrants are weighted by the inverse of their quantity.
 - 1/X^2: Specifies a weighting of 1/X^2 for all calibration data points during the least-squares regression calculation of the calibration curve. Calibrants are weighted by the inverse of the square of their quantity.
 - 1/Y: Specifies a weighting of 1/Y for all calibration data points during the least-squares regression calculation of the calibration curve. Calibrants are weighted by the inverse of their response (or response ratio).
 - 1/Y^2: Specifies a weighting of 1/Y^2 for all calibration data points during the least-squares regression calculation of the calibration curve. Calibrants are weighted by the inverse of the square of their response (or response ratio).
- 5. From the Response Via list, select Area or Height.
 - Area: Specifies that the TraceFinder application use this area value in response calculations.
 - Height: Specifies that the application use this height value in response calculations.

To specify qualitative peak processing



1. Select the **Use Genesis Algorithm for Qual Processing** check box and specify a value for internal standard matching.

The application uses the Genesis algorithm to match internal standards in a range plus/minus the value you specify. For additional information about the Genesis algorithm, see "Genesis Detection Method" on page 73.

2. Select or clear the **Exclude Matching Quan Peaks** check box and specify a value for the exclusion window.

The application excludes quan peaks in a range plus/minus the value you specify.

3. To process samples that include data-dependent scans, select the **Use Data Dependent Scans** check box.

When you process a sample using this feature, the application uses the TIC trace to find all data-dependent full scans, lists them, and performs a library search against the data-dependent MS/MS or MSⁿ scan.

In addition to the peak information, the TIC Report and TIC Summary Report display information about the data-dependent filtered data. See Appendix A, "Reports."

Working with Master Methods

To enter a note for the method

Type in the Notes box, or paste text from another application using CTRL+V.

You can add a note to your method template to explain what makes this template unique.

❖ To save the method template

1. Choose **File > Save** from the Method Template Editor menu.

The Save Method Template dialog box opens.

2. Do one of the following:

Type a new name for the master method and click **OK**.

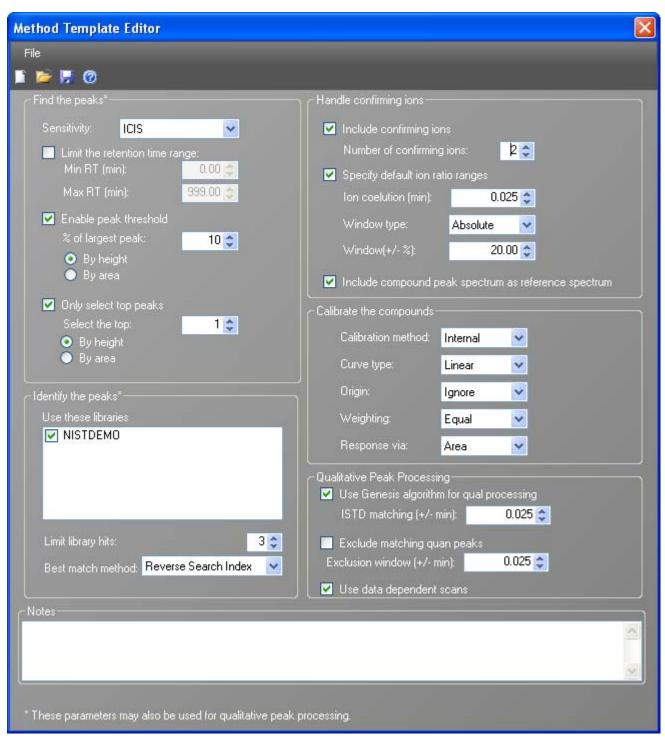
-Or-

Select a method name to overwrite and click **Overwrite**.

The TraceFinder application saves the new method template in the following folder:

 $... \label{thm:linear} TraceFinder \verb|\| 1.1 \\ Templates \\ Methods$

Figure 58. Method Template Editor dialog box



Working with Master Methods

Table 52. Method Template Editor dialog box parameters (Sheet 1 of 3)

| Parameter | Description |
|--|---|
| Find the peaks | |
| Sensitivity | Defines how extensively the peak detector algorithm searches for low-level peaks. |
| Limit the retention time range | Min RT specifies the beginning of the range. Max RT specifies the end of the range. |
| Enable peak threshold | Specifies whether to select peaks by relative height or area and the percentage of the highest peak that results in compound selection. |
| Only select top peaks | Displays a specific number of the largest peaks by height or area. |
| Identify the peaks | |
| Use these libraries | Lists the libraries you can search. |
| Limit library hits | Specifies the number of hits returned when the system searches a spectrum against the selected libraries. |
| Best match method | Specifies how to sort the library searches. Valid values: Search Index, Reverse Search Index, Match Probability |
| Handle confirming ions | |
| Include confirming ions/ Number of confirming | Specifies the number of confirming ions, which are other ions in the spectrum whose ratio is compared to the quantitation ion to identify the compound. |
| ions | This value defaults to 2 because you typically perform a 3-ion experiment with one quan mass and two confirming ions. |
| Specify default ion ratio | Enables the ion ratio range features. |
| ranges | Ion Coelution specifies the maximum difference in retention time between a confirming ion peak and the quantification ion peak. |
| | Window Type specifies an Absolute or Relative calculation approach for determining the acceptable ion ratio range. |
| | Window (+/-%) specifies the acceptable ion ratio range. |
| Include compound peak spectrum as reference spectrum | Includes the peak spectrum in the processing method. Use this setting to perform a spectra comparison in Data Review. |
| Calibrate the compounds | |
| Calibration method | Specifies an internal or external calibration method. |
| Curve type | Specifies a linear, quadratic, or average RF curve type. |
| * 1 | |

Table 52. Method Template Editor dialog box parameters (Sheet 2 of 3)

| Parameter | Description |
|---|---|
| Origin | Specifies that the origin is ignored, forced, or included in the generated calibration curve. Ignore: Specifies that the origin is not included as a valid point in the calibration curve when the curve is generated. When you select Ignore, the calibration curve might or might not pass through the origin. Force: Specifies that the calibration curve passes through the origin of the data point plot when the calibration curve is generated. Include: Specifies that the origin is included as a single data point in the calculation or the calibration curve. When you select Include, the calibration curve might or might not pass through the origin. |
| Weighting | Specifies the weighting for the calibration data points. Equal: Specifies that the origin is included as a single data point in the calculation of the calibration curve. When you select Equal, the calibration curve might or might no pass through the origin. 1/X: Specifies a weighting of 1/X for all calibration data points during the least-squares regression calculation of the calibration curve. Calibrants are weighted by the inverse of their quantity. 1/X^2: Specifies a weighting of 1/X^2 for all calibration data points during the least-squares regression calculation of the calibration curve. Calibrants are weighted by the inverse of the square of their quantity. 1/Y: Specifies a weighting of 1/Y for all calibration data points during the least-squares regression calculation of the calibration curve. Calibrants are weighted by the inverse of their response (or response ratio). 1/Y^2: Specifies a weighting of 1/Y^2 for all calibration data points during the least-squares regression calculation of the calibration curve. Calibrants are weighted by the inverse of their response of their response ratio). |
| Response via | Specifies if the TraceFinder application uses area or height in response calculations. Area: Specifies that the application use this peak area value in response calculations. Height: Specifies that the application use this peak height value in response calculations. |
| Qualitative Peak Process | ing |
| Use Genesis algorithm for qual processing | The application uses the Genesis algorithm to match internal standards. |
| ISTD matching | Excludes all the target compounds found in the method and does not list these compounds in the TIC Report or in the Qual Mode view in the Data Review. |

Table 52. Method Template Editor dialog box parameters (Sheet 3 of 3)

| Parameter | Description |
|-----------------------------|---|
| Exclude matching quan peaks | Compares the retention time of the internal standard in the method to the found retention time of the internal standard in the library search and excludes peaks outside the Exclusion Window range. |
| Exclusion window | Defines a range plus/minus the Exclusion Window value you specify. |
| Use data dependent scans | Constrains the Qual Mode view in the Data Review to only data-dependent scan spectra. See "Qual Mode" on page 313. In addition to the peak information, the TIC Report and TIC Summary report display information about the data-dependent filtered data. |

Importing Published Master Methods

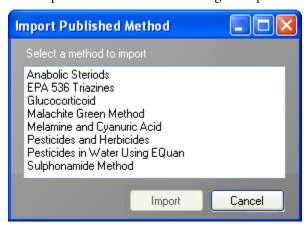
In the TraceFinder application, you can import published methods to use for detecting, processing, and reporting. The Tracefinder installation provides the following folder of published methods:

...\Thermo\TraceFinder\1.1\Published Master Methods

❖ To import a published master method

1. From the Method View task pane, click **Import Published Method**.

The Import Published Method dialog box opens.



2. Select a method to import.

3. Click **Import**.

The application reports that the method successfully imported and saves the method in the following folder:

 $...\Thermo\TraceFinder\1.1\Methods$

You can use any of the Open Method commands to open this method just as you would a method you created.

Exporting SRM Data

In the TraceFinder application, you can export your selected reaction monitoring (SRM) data to an XML file. The Export SRM Data command is available only when you select the Enable Compound Datastore option on the General page in the Application Configuration view. See "Application Configuration" on page 56.

❖ To export SRM data to an XML file

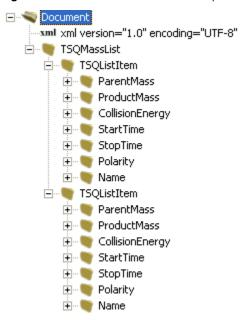
- 1. Open the master method whose SRM data you want to export.
- 2. From the Method View task pane, click Export SRM Data.

The TraceFinder application writes the data in the SRM table to the following file:

...\Thermo\TraceFinder\1.1\Methods\methodname.xml

The data in this file matches the TSQ .xml data, which you can use in the instrument method editor of the TSQ application.

Figure 59. SRM TSQ Quantum™ example



Working with Instrument Methods

An instrument method is a set of experiment parameters that define the operating settings for an autosampler, mass spectrometer, and so on. Instrument methods are saved as file type .meth.

IMPORTANT Do not open the Thermo Foundation Instrument Configuration window while the TraceFinder application is running.

Use the following procedures:

- To open the Instrument View
- To create a new instrument method
- To create a new multiplexing instrument method
- To open an instrument method
- To import an instrument method
- To open the Instrument View
- 1. Click **Method Development** from the dashboard or the navigation pane.



The Method Development navigation pane opens.

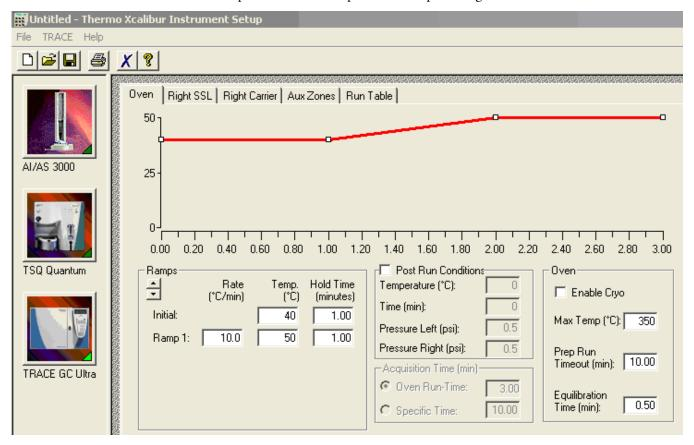
2. Click the **Instrument View** task pane.



1. Click New Instrument Method in the Instrument View task pane.

The Thermo Xcalibur Instrument Setup window opens.

This example instrument setup shows multiple configured instruments.



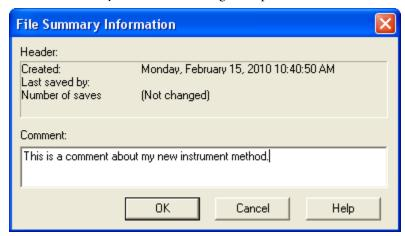
- 2. Click the icon for the instrument you want to use for the method.
- 3. Edit the values on the instrument page.
- 4. From the main menu in the Thermo Xcalibur Instrument Setup window, choose **File > Save As**.

The Save As dialog box opens.

Working with Instrument Methods

5. Select an instrument method name to overwrite or type a new name for the instrument method, and click **Save**.

The File Summary Information dialog box opens.



- 6. (Optional) Type a comment about the new instrument method.
- 7. Click OK.

The TraceFinder application saves the new instrument method in the following folder:

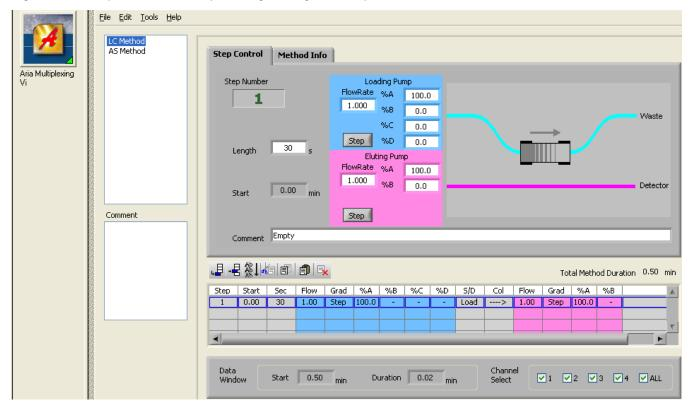
...\Xcalibur\methods

❖ To create a new multiplexing instrument method

1. Click New Instrument Method in the Instrument View task pane.

The Thermo Xcalibur Instrument Setup window opens.

Figure 60. Example instrument setup showing a configured multiplexed instrument



- 2. Click the icon for the instrument you want to use for the method.
- Edit the values for the instrument method.
 For information about specifying multiplexing values, refer to the documentation for your multiplexed instrument.
- 4. Specify the channels you want to use for acquisition. For example:



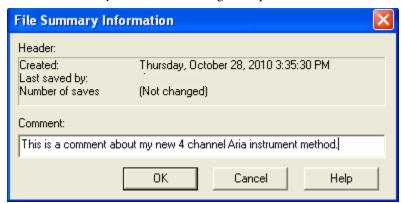
5. From the main menu in Thermo Xcalibur Instrument Setup window, choose **File > Save As**.

The Save As dialog box opens.

Working with Instrument Methods

6. Select an instrument method name to overwrite or type a new name for the instrument method, and click **Save**.

The File Summary Information dialog box opens.



- 7. (Optional) Type a comment about the new instrument method.
- 8. Click OK.

The TraceFinder application saves the new instrument method in the following folder:

 $...\X$ calibur\methods

❖ To open an instrument method

1. Click **Open Instrument Method** on the Instrument View task pane.

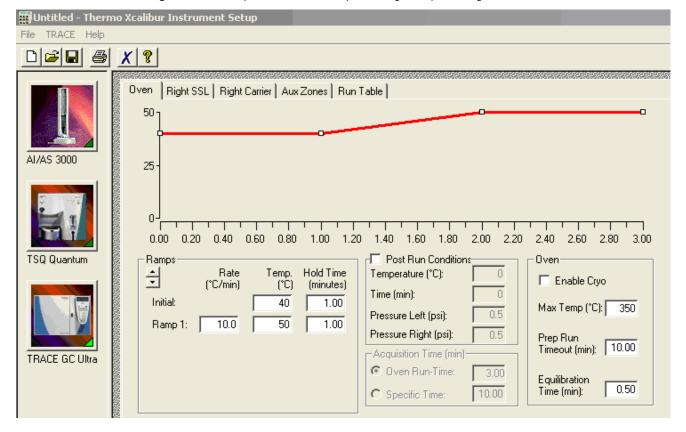
An instrument method browser opens.

2. In the browser, select an instrument method from the list and open the file.

The selected method opens in the Thermo Xcalibur Instrument Setup window. You can edit this method and save the changes, or you can save this method to another name.

Note To open Help for any of your configured instruments, click **Help** on the instrument page.

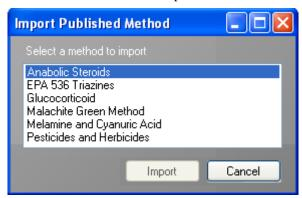
Figure 61. Example instrument setup showing multiple configured instruments



❖ To import an instrument method

1. From the Instrument View task pane, click **Import Published Method**.

The Import Published Method dialog box opens. This dialog box lists the master methods in the Published Master Methods folder. You can import instrument methods that are associated with these published master methods.

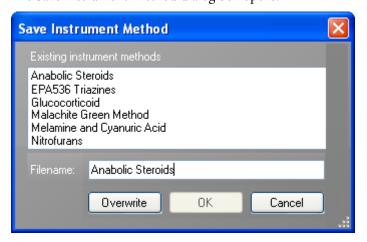


2. Select a method that includes the instrument method you want to import.

For instructions for importing the master methods, see "Importing Published Master Methods" on page 198.

3. Click **Import**.

The Save Instrument Method dialog box opens.



4. Do one of the following:

Type a new name for the instrument method and click **OK**.

-Or-

Select an instrument method name to overwrite and click **Overwrite**.

The application reports that the method successfully imported.

You can use any of the Open Instrument Method commands to open this method just as you would an instrument method that you created.

Working with Development Batches

In the Development Batch view, you can test your instrument method in real time by creating and acquiring test samples. Development batches let you test different instrument methods and optimize parameters, such as MS source parameters and autosampler variables, to find the best conditions for a master method. Development batches are not designed for high throughput in everyday analysis.

This section includes instructions for the following tasks:

- Creating a Development Batch
- Editing Samples in a Development Batch
- Acquiring Samples in a Development Batch

Creating a Development Batch

You create a development batch to test your instrument method and use it to acquire samples only once. You cannot save a development batch; you can save only the raw data files created when you acquire the samples in the batch.

Use the following procedures:

- To open the Development Batch view
- To specify a location for development batch data
- To add samples to the development batch
- To insert samples into the development batch
- To copy a sample

To open the Development Batch view

1. Click **Method Development** from the dashboard or the navigation pane.



The Method Development navigation pane opens.

2. In the Method Development navigation pane, click **Development Batch**.



The Development Batch view opens a new, empty batch.



Note The Channel column is available only when your configuration supports multiplexing.

To specify a location for development batch data

1. To specify a location for the files, click **Select Batch Location** in the Development Batch task pane.

By default, the TraceFinder application writes the temporary files, raw data files, and .sld method file to the following folder:

- ...\Thermo\TraceFinder\1.1\Temp
- 2. In the browser, do one of the following:

Locate the folder you want to use for the development batch files and click **OK**.

-Or-

Do the following:

- a. Locate and select the folder where you want to create a new folder for the batch files
- b. Click Make New Folder.

The TraceFinder application creates a new folder in the selected folder.

- c. Right-click the New Folder filename and choose **Rename** from the shortcut menu.
- d. Type the name for the folder.
- e. Click **OK**.

The TraceFinder application creates all development batch files in the specified folder.

❖ To add samples to the development batch

Do one of the following:

Right-click and choose **Add Sample** from the shortcut menu.

-Or-

To add multiple sample rows, enter the number of rows and 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 development batch

- 1. Select the sample above which you want to insert empty samples.
- 2. Do one of the following:

Right-click and choose Insert Sample from the shortcut menu.

-Or-

To insert multiple sample rows, enter the number of rows and click the **Insert Sample** icon.



The TraceFinder application inserts new, empty samples above the selected sample.

Note You cannot insert samples into an empty batch. You must have at least one sample to select before you can use this icon.

❖ To copy a sample

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

The TraceFinder application adds a copy of the sample above the selected sample.

Editing Samples in a Development Batch

A development batch requires fewer parameters than a real batch, but the mechanism for managing the information is the same.

For detailed instructions about using the Copy Down or Fill Down commands to enter column values, see Appendix B, "Using Copy Down and Fill Down."

Use the following procedures:

- To enter column values
- To resize or reorganize the columns
- To remove selected samples from the list
- To remove all samples from the list

❖ To enter column values

- 1. Double-click the Filename column and type a file name for the raw data file.
- 2. (Optional) Enter values for the Sample Name or Sample ID columns.
- 3. Enter a vial position for each sample.
- 4. Enter an injection volume for each sample.

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

5. To enter an instrument method for each sample, click the down arrow in the Instrument Method column and select a method from the list.

This list contains all the available instrument methods.

6. To enter a channel for each sample, click the down arrow in the Channel column and select a channel from the list.



You cannot specify the auto channel selection in a development batch.

Note The Channel column is available only when your configuration supports multiplexing.

Figure 62. Completed Development Batch

| Development Batch - [C:\Thermo\TraceFinder\1.1\Temp]* | | | | | | | | |
|---|---|-------|---|--|---|------|---------------------|-----------|
| Filename Sample Sample Vial Injection Instrument Channe | | | | | | | | |
| | 1 | File1 | 1 | | 1 | 10.0 | Anabolic Steroids 🔽 | Channel 1 |
| | 2 | File2 | 2 | | 2 | 10.0 | Anabolic Steroids 🔽 | Channel 2 |
| | 3 | File3 | 3 | | 3 | 10.0 | Anabolic Steroids 🔽 | Channel 1 |
| • | 4 | File4 | 4 | | 4 | 10.0 | Anabolic Steroids 🔽 | Channel 2 |

❖ To resize or reorganize the columns

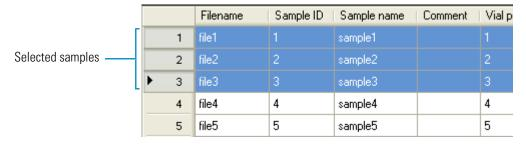
- 1. To resize a column, drag the header separator on the right side of the column.
- 2. To move a column, drag the column header.

You cannot move the Filename column.

❖ To remove selected samples from the list

1. Select the samples you want to remove.

Use the first column to ensure that the samples are selected.



2. Right-click and choose **Remove Selected Samples** from the shortcut menu.

❖ To remove all samples from the list

1. Click New Sample List in the Development Batch task pane.

One of the following happens:

- If the samples in the current batch have all been acquired, the list is cleared.
- If the samples in the current list have not been acquired, a message confirms that you want to clear them and start a new list.
- 2. To create a new empty list, click Yes.

Note You cannot save a development batch when you create a new one; you can only create, acquire, and discard each batch after you use it. The TraceFinder application saves only the generated raw data files in the specified batch location.

Acquiring Samples in a Development Batch

In a development batch, you can submit the entire batch for acquisition or submit only selected samples.

Use the following procedures:

- To acquire selected samples
- To acquire the batch

❖ To acquire selected samples

- 1. Select the samples you want to acquire.
- 2. Right-click and choose **Submit Selected Samples** from the shortcut menu, or click the

Submit Selected Samples icon,

The TraceFinder application creates a raw data file for each selected sample. It writes the raw data files and all temporary working files to the following folder:

...\Thermo\TraceFinder\1.1\Temp

When the acquisition is complete, the application deletes all the temporary working files. Only the raw data files and a MethodDevelopment.sld file remain in the folder.

If a sample is acquired more than once, the subsequent raw data files are time-stamped with the acquisition time.

To acquire the batch

Right-click and choose **Submit Batch** from the shortcut menu, or click the **Submit Batch** icon,

The TraceFinder application creates a raw data file for each sample in the batch and an .sld method file. The TraceFinder application writes the raw data files, the .sld method file, and all temporary working files to the specified folder.

When the acquisition is complete, the application deletes all the temporary working files. Only the raw data files and a MethodDevelopment.sld file remain in the folder.

If a sample is acquired more than once, the subsequent raw data files are time-stamped with the acquisition time.

Viewing Raw Data Files in the Qual Browser

You can view the chromatogram and spectra for completed samples in a development batch.

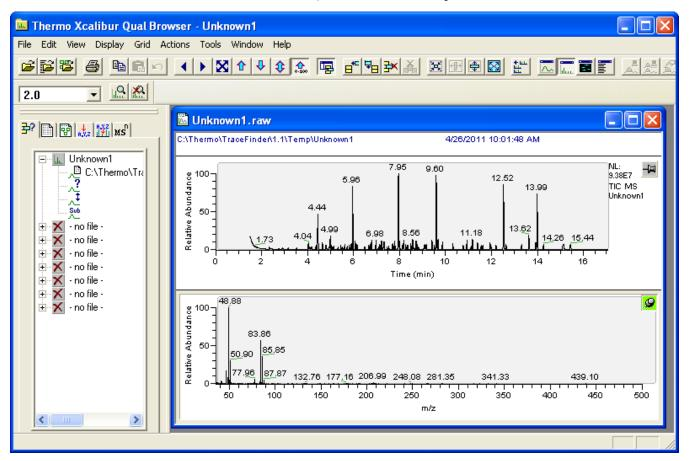
Use the following procedures:

- To open the Qual Browser
- To display the last completed raw data file in the Qual Browser

❖ To open the Qual Browser

In the Development Batch task pane, click Open Qual Browser.

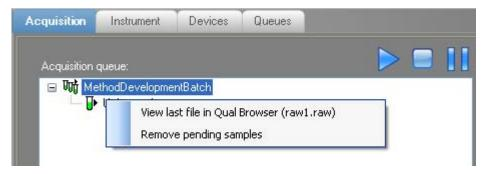
The Thermo Xcalibur Qual Browser window opens.



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

❖ To display the last completed raw data file in the Qual Browser

In Acquisition page of the real-time viewer, right-click and choose **View Last File in Qual Browser** from the shortcut menu.



The last completed file opens in the Qual Browser.

When all samples are completed, you can view the last raw data file for the batch.



Using the Acquisition Mode

This chapter describes the tasks associated with the Acquisition mode.

Contents

- Working with Batches
- Real-time Display
- 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), a user in the LabDirector or Supervisor role can define a batch template that supplies the basic structure of a batch.

If you have a master method, you can create a batch and run the samples. Batches represent 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 the Data Review mode or you can go directly to viewing and printing reports.

You can also set up a calibration batch with known concentrations of the target compounds to compare the sample against the calibration in future batches.

Working with Batches

This section includes instructions for the following tasks:

- Opening and Navigating the Acquisition Mode
- Creating Batches

Opening and Navigating the Acquisition Mode

To open the Acquisition mode, click **Acquisition** in the navigation pane or from the dashboard.

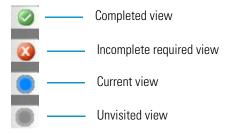


The TraceFinder application does not use the navigation pane in the Acquisition mode in the same way it uses the navigation pane in other modes. In the Acquisition mode, this pane keeps track of your progress as you move through the views to create and submit a batch or a batch template.

Figure 63. Task pane when you enter the Acquisition mode

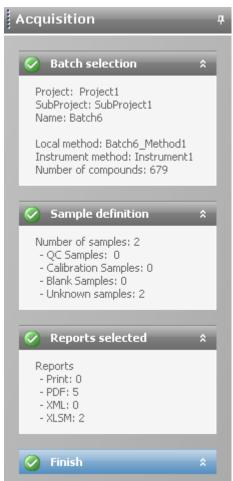


The status of each view in the Acquisition mode shows you which tasks are completed and which tasks are not.



As you complete each view, the task panes display the parameters you specified for your batch.

Figure 64. Example task pane when you have completed the Acquisition mode



Creating Batches

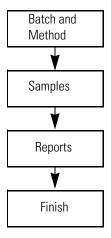
These are the major steps you use in Acquisition mode to create a batch:

- 1. Selecting a Batch
- 2. Defining the Sample List
- 3. Selecting and Reviewing Reports
- 4. Submitting the Batch

Workflows

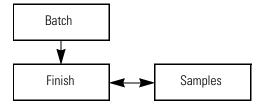
Depending on how you create your batch, you use different views in the Acquisition mode. The following workflows show the Acquisition views required for each batch creation approach. Depending on your approach to creating a batch, use one of these specific workflows.

Creating an original batch



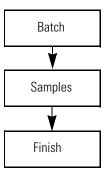
To create an original batch, start with the instructions "To start a new batch" on page 220.

❖ Acquiring a previously saved (.tbr) batch



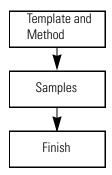
To acquire a previously saved batch, start with the instructions "To select a ready-to-acquire batch" on page 221.

& Editing and processing a previously acquired batch



To process a previously acquired batch, start with the instructions "To select a previously acquired batch" on page 222.

❖ Creating a batch template



When user security is enabled, this workflow is available only to users in the LabDirector or Supervisor role.

To create a batch template, start with the instructions "To create a batch template" on page 223 and then click **Save** on the Finish page.

Selecting a Batch

In a Template Selection view of the Acquisition mode, you can choose to create a new batch in any of your current projects/subprojects.

Use the following procedures:

- To start a new batch
- To select a template for the new batch
- To select a ready-to-acquire batch
- To select a previously acquired batch
- To create a batch template

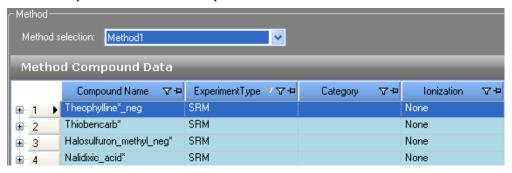
❖ To start a new batch

- 1. Click the **New Batch** tab.
- 2. Select the project and subproject where you want to create the new batch.
- 3. Type a name for the new batch in the Batch Name box.



4. 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.



5. To continue to the next view, click **Next**.

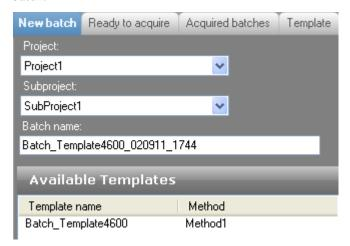
The Sample Definition view opens. See "Defining the Sample List" on page 224.

To select a template for the new batch

- 1. Click the New Batch tab.
- 2. In the Available Templates pane, select the template and method combination you want to use.

The system creates a Batch name with the selected template name and the date/time stamp. You can change the default project, subproject, and method associated with this template.

3. (Optional) Select a different project and subproject where you want to create the new batch.



- 4. (Optional) Select a different method to use for the new batch.
- 5. To continue to the next view, click **Next**.

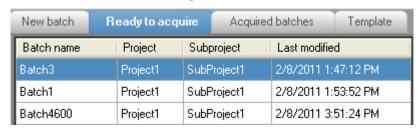
The Sample Definition view of the Acquisition mode opens. See "Defining the Sample List" on page 224.

To select a ready-to-acquire batch

1. Click the **Ready to Acquire** tab.

All your unacquired, saved batches are displayed with the file extension .tbr (to be run). All .tbr batches are located in the following folder:

- ...\Thermo\Tracefinder\1.1\Projects\projectname\subprojectname
- 2. Select the batch you want to acquire.



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

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

- 4. Do one of the following:
 - To prepare the batch for acquisition, click **Submit**.
 For detailed instructions, see "Submitting the Batch" on page 237.
 - -Or-
 - To edit the samples list, click **Previous**.

For detailed instructions on editing a samples list, see "Defining the Sample List" on page 224.

-Or-

• To save the batch to the Ready to Acquire list, click **Save**.

The TraceFinder application saves your batch as a to-be-run (.tbr) file, closes the Acquisition mode, and returns you to the application dashboard.

The application saves the .tbr batches in the following folder:

...\Thermo\TraceFinder\1.1\Projects\projectname\subprojectname

To select a previously acquired batch

1. Click the Acquired Batches tab.

From this page, you can resubmit a previously acquired batch, edit the batch, or save it to be acquired later.

2. In the Project pane, select a project name.

All subprojects included in the selected project are displayed in the Subproject pane.

3. In the Subproject pane, select a subproject name.

All previously acquired batches included in the selected subproject are displayed in the Batch pane.

4. In the Batch pane, select the batch you want to reacquire.



5. To continue to the next view, click **Next**.

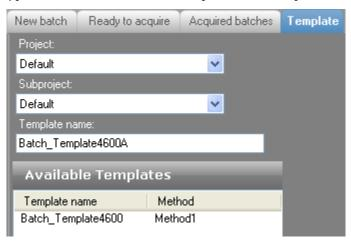
The Sample Definition view of the Acquisition mode opens. See "Defining the Sample List" on page 224.

To create a batch template

1. Click the **Template** tab.

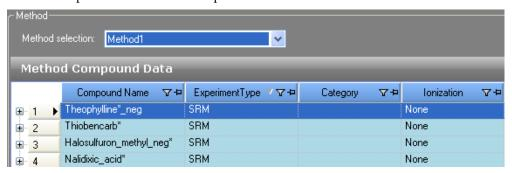
Note When user security is enabled, this page is available only to users in the LabDirector or Supervisor role.

- 2. Select the project and subproject where you want to create the new batch template.
- 3. Type a name for the new batch template in the Template Name box.



4. 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.



5. To continue to the next view, click **Next**.

The Sample Definition view of the Acquisition mode opens. See "Defining the Sample List" on page 224.

Defining the Sample List

In the Sample Definition view of the Acquisition mode, you can create a list of samples for the batch. You can add samples, insert samples, import a sample list, or remove samples from the list. See "Sample Definition view" on page 230.

As you enter sample values, you can use the Copy Down and Fill Down commands to 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 the following 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

When you have finished defining the list of samples, click Next.

- When you are creating a batch from scratch, creating a batch from a template, or editing
 a batch template, the Report Selection view opens. See "Selecting and Reviewing Reports"
 on page 233.
- When you are editing a previously acquired batch or a .tbr batch, the Finish Selection view opens. See "Submitting the Batch" on page 237.

To add samples to the list

1. Select the number of sample rows to add and click the **Add** button, Add



Tip To quickly add a single row, right-click and choose **Add Sample** from the shortcut menu.

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 Environmental and Food Safety sample types | | | | | |
|--|------------|-------------|--|--|--|
| Matrix Blank | Solvent | Unknown/TIC | | | |
| Cal Std | Chk Std | Unknown | | | |
| LCS | MDL | MS | | | |
| LCSD | Method Val | MSD | | | |

| Available General Quantitation sample types | | | | |
|---|-------------|---------|--|--|
| Matrix Blank | Unknown/TIC | Unknown | | |
| Cal Std | Chk Std | Solvent | | |

For a detailed description of sample types, see "Sample Types" on page 259.

4. For each Cal Std or Chk Std sample, select a level from the Sample Level list.

The sample levels are defined in the master method. If there are no levels to select from the Sample Level list, ask a user with Supervisor or LabDirector permissions to edit the method and specify the levels.

For detailed instructions about defining sample levels, see Chapter 4, "Using the Method Development Mode."

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

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

6. 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 .

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

Note When you use the horizontal scroll bar at the bottom of the samples list, the Status and Filename columns stay fixed while the other columns scroll right and left.

❖ To insert samples into the list

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

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



The application inserts the unknown samples above the selected sample.

| | | Status | Filename | Sample type | Sample level | Sample ID |
|--------------------|---|--------|------------|-------------|--------------|--------------------|
| Inserted samples – | 1 | | cal_std_5 | Cal Std | 10 | cal std = 5 ng/uL |
| | 2 | | Unknown2 | Unknown | | |
| | 3 | • | Unknown1 | Unknown | | |
| | 4 | • | cal_std_10 | Cal Std | 10 | cal std = 10 ng/uL |

3. Type a file name in the Filename column for each sample.

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

Each file name must be unique.

4. Select a sample type from the Sample Type list for each sample.

For a detailed description of each sample type, see "Sample Types" on page 259.

| Available Environmental and Food Safety sample types | | | | |
|--|------------|-------------|--|--|
| Matrix Blank | Solvent | Unknown/TIC | | |
| Cal Std | Chk Std | Unknown | | |
| LCS | MDL | MS | | |
| LCSD | Method Val | MSD | | |

| Available General Quantitation sample types | | | | | |
|---|-------------|---------|--|--|--|
| Matrix Blank | Unknown/TIC | Unknown | | | |
| Cal Std | Chk Std | Solvent | | | |

5. For each Cal Std or Chk Std sample, click the Sample Level cell and select a level from the list.

The sample levels are defined in the master method. If there are no levels to select from the Sample Level list, ask a user with Supervisor or LabDirector permissions to edit the method and specify the levels.

For detailed instructions about defining sample levels, see Chapter 4, "Using the Method Development Mode."

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 μ L; the maximum injection volume value allowed is 5000 μ L.

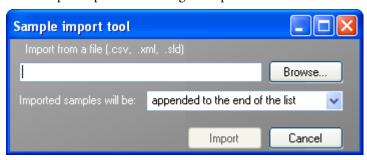
8. (Optional) Type or edit the values for the remaining columns.

Note When you use the horizontal scroll bar at the bottom of the samples list, the Status, Filename, Sample Type, and Sample Level columns stay fixed while the other columns scroll right and left.

❖ To import samples into the list

1. Click **Import**, Import

The Sample Import Tool dialog box opens.



From this dialog box, you can import a sample list from a .csv, .xml, or .sld file.

2. Click **Browse** and select a .csv, .xml, or .sld file that contains the sample definitions you want to import.

Note The .csv, .xml, or .sld file format must match the TraceFinder file 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 specified samples are added to the Samples list.

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

| Xcalibur column | TraceFinder column |
|-----------------|--------------------|
| Level | Sample Level |
| Position | Vial Position |
| Inj Vol | Injection Volume |
| Dil Factor | Conv Factor |

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

| Xcalibur sample type | TraceFinder sample type |
|----------------------|-------------------------|
| Blank | Matrix Blank |
| QC | Chk Std |
| Std Bracket | Cal Std |

5 Using the Acquisition Mode

Working with Batches

5. For each Cal Std or Chk Std sample, click the Sample Level cell and select a level from the list.

The sample levels are defined in the master method. If there are no levels to select from the Sample Level list, ask a user with Supervisor or LabDirector permissions to edit the method and specify the levels.

For detailed instructions about defining sample levels, see Chapter 4, "Using the Method Development Mode."

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 horizontal scroll bar at the bottom of the samples list, the Status, Filename, Sample Type, and Sample Level columns stay fixed while the other columns scroll right and left.

(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 you want to remove.

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

2. Right-click and choose **Remove Selected Samples** from the shortcut menu.

To reinject a sample from a previously acquired batch

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

The TraceFinder 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 TraceFinder 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).



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

❖ To select channels for the batch

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



By default, all configured channels are selected. The configured channels are determined by the multiplexing settings in the Configuration mode. See "Specifying Multiplexing Parameters" on page 70.

Disabling 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 you disabled.

❖ To assign a specific channel to a sample

- Scroll to the Channel column (the rightmost column in the sample list).
 All samples default to Auto.
- 2. Select a channel from the Channel list.

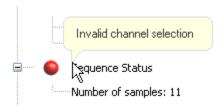


When you submit the batch, samples that are set to Auto run on any of the available channels and samples that are set to a specific channel run only on that channel.

5 Using the Acquisition Mode

Working with Batches

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.

Figure 65. Sample Definition view



Table 53. Sample Definition parameters (Sheet 1 of 3)

| Parameter Definition | | | |
|-----------------------|---|--|--|
| 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 where you can import samples defined in a .csv file or an .xml file. | | |
| Multiplexing Channels | | | |
| All Channels | Uses all configured channels to acquire this batch. | | |
| Channel 1-n | Uses only the selected channels to acquire this batch. | | |
| Previous | Returns you to the previous Acquisition mode view. | | |
| Cancel | Confirms that you want to exit the Acquisition mode. When you cancel out of the Acquisition mode, your edits are not saved. | | |
| Save | Saves this batch as a to-be-run (.tbr) batch. | | |

Table 53. Sample Definition parameters (Sheet 2 of 3)

| Parameter | Definition | | | | |
|---------------------------|---|--|--|--|--|
| Next | Takes you to the next Acquisition mode view. | | | | |
| Shortcut menu | | | | | |
| 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" on page 227. | | | | |
| 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" on page 286. | | | | |
| Сору | Copies the data in the selected rows or columns to the Clipboard. Use this command to copy sample information to another application, such as an Excel spreadsheet. 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 another application, such as an Excel spreadsheet. You cannot paste this data back into the Acquisition mode sample list. For example: | | | | |
| | Sample type Matrix Blank Cal Std Chk Std Unknown/TIC Unknown/TIC Unknown/TIC Copy With Headers from TraceFinder | | | | |
| Paste | Pastes a single column of copied data from another application, such as an Excel spreadsheet, into the selected column. | | | | |
| Undo Last Paste | Removes the last pasted item in the Acquisition mode sample list. | | | | |

5 Using the Acquisition Mode Working with Batches

Table 53. Sample Definition parameters (Sheet 3 of 3)

| Parameter | Definition |
|--------------------|---|
| Export to CSV file | Opens the Save As dialog box where you can save the current sample list to a .csv file. |
| Status color codes | Sample is not acquired. |
| | Sample is acquired but not processed. |
| | Sample is acquired and processed. |

Selecting and Reviewing Reports

On the Report Selection view, you can specify the types of reports you want to create. See "Report Selection view" on page 235. For a complete list of report types and examples of output files, see Appendix A, "Reports." In addition to the report type, you can specify a report description for each of your reports.

For each standard report you generate, you can create a hardcopy printout, a PDF file, or an XML file.

For each custom report you generate, you can create a hardcopy printout or an XLSM file.

When you have finished specifying your report options, click **Next** to go to the Finish view. See "Submitting the Batch" on page 237.

The resulting output files for your reports are written to the following folder:

...\Projects\projectname\subprojectname\batchname\Reports

Use the following procedures:

- To edit a report description
- To preview a standard report
- To specify a standard report in print format or as a PDF, XML, or XLSM file
- To export reports to a specific folder

❖ To edit a report description

Select the Report Description column and edit the default description.

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

❖ To preview a standard report

Click the magnifying icon, , to view an example of the report type as a PDF file.
 The right pane of the view displays an example PDF report with typical PDF viewer buttons.

2. To minimize the PDF viewer, click the minimize icon,

Note Only Standard report types have preview documents.

* To specify a standard report in print format or as a PDF, XML, or XLSM file

- 1. For each type of report you want to create, select the check box in the Print, Create PDF, Create XML, or Create XLSM columns.
- 2. To duplicate the output type for all reports, right-click the cell and choose **Copy Down** from the shortcut menu.

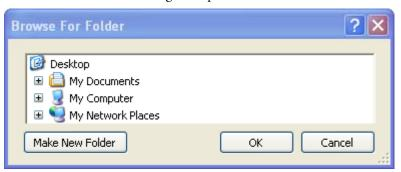
All check boxes in the column below the selected cell duplicate the selected or cleared state in the selected cell. You can duplicate the output type only for reports that have this output format available.

❖ To export reports to a specific folder

1. Select the **Export Results** check box at the bottom of the view.



The Browse For Folder dialog box opens.



- 2. Locate and select the folder where you want to save the reports.
- 3. To create a new reports folder within the selected folder, click **Make New Folder** and type the new folder name.
- 4. Click OK.

The application writes all reports to the specified folder in addition to the batch Reports folder.

Figure 66. Report Selection view

| Example | Report Name | Report Description | Report Type | Print | Create PDF | Create XML | Create XLSM | Batch Lev |
|---------|-----------------------------|-----------------------------|----------------|-------|------------|------------|-------------|-----------|
| ۶ | Batch Report | Batch Report | Standard | | | | | |
| P | Blank Report | Blank Report | Standard | | | | | |
| P | Calibration Report | Calibration Report | Standard | | | | | |
| P | Chromatogram Report | Chromatogram Report | Standard | | | | | |
| P | Compound Calibration Re | Compound Calibration Re | Standard | | | | | |
| P | Compound Calibration Re | Compound Calibration Re | Standard | | | | | |
| P | Confirmation Report | Confirmation Report | Standard | | | | | |
| P | Confirmation Report 2 | Confirmation Report 2 | Standard | | | | | |
| P | High Density Calibration R | High Density Calibration R | Standard | | | | | |
| P | High Density Internal Stan | High Density Internal Stan | Standard | | | | | |
| P | High Density Internal Stan | High Density Internal Stan | Standard | | | | | |
| P | High Density Sample Rep | High Density Sample Rep | Standard | | | | | |
| P | High Density Sample Rep | High Density Sample Rep | Standard | | | | | |
| P | High Density Sample Rep | High Density Sample Rep | Standard | | | | | |
| P | High Density Sample Rep | High Density Sample Rep | Standard | | | | | |
| P | AltCalibrationReport | AltCalibrationReport | Custom | | | | | |
| P | Alternate BatchReport | Alternate BatchReport | Custom | | | | | |
| P | Alternate CalibrationReport | Alternate CalibrationReport | Custom | | | | | |
| P | Alternate ConfirmationRep | Alternate ConfirmationRep | Custom | | | | | |
| P | Alternate MatrixSpikeReport | Alternate MatrixSpikeReport | Custom | | | | | |
| P | Alternate SampleReport | Alternate SampleReport | Custom | | | | | |
| P | Alternate SummaryReport | Alternate SummaryReport | Custom | | | | | |

Table 54. Report Selection parameters (Sheet 1 of 2)

| Parameter | Description | |
|--------------------|--|--|
| P | Displays an example PDF for the report type. This example provides a model of the report type only; it does not reflect your specific data. This is available for standard reports only. | |
| Report Name | The name of a report. | |
| Report Description | User-editable title to be used on a report. | |
| Report Type | The type of report: Standard, Custom, or Target Screening. | |
| Print | Reports to be sent to the printer. | |
| Create PDF | Reports to be saved as PDF files. Available only for standard reports. | |
| Create XML | Reports to be exported in XML format. Available only for standard reports. | |

5 Using the Acquisition Mode Working with Batches

 Table 54. Report Selection parameters (Sheet 2 of 2)

| Parameter | Description |
|-----------------------------|--|
| Create XLSM | Reports to be exported in Excel Macro-Enabled Workbook (.xlsm) format. Available only for custom reports. |
| Batch Level | Rather than creating separate reports for each sample, the application uses a composite of the data from all the samples to create a single report for the entire batch. Batch-level reports are prepended with a B to differentiate them. You cannot select this option from the Report Selection page. You must select the Batch Level option for the report in the report configuration. See "Specifying the Reports Configuration" on page 58. |
| Shortcut menu: Copy Down | Copies the selected or cleared state to all subsequent reports in the column. |

Submitting the Batch

In the Finish view 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. See "Finish view" on page 244.

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

Use the following procedures:

- To specify startup or shutdown methods
- To specify a calibration batch
- To specify device states
- To save a batch for later acquisition
- To start an acquisition
- To view the output files

To specify startup or shutdown methods

1. (Optional) Select a method from the System Startup Method list.

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

2. (Optional) Select a method from the System Shutdown Method list.

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



5 Using the Acquisition Mode

Working with Batches

❖ 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 .calx calibration file.

2. Do one of the following:

To use the selected calibration file to process the current data, select the **Use Calibration** option.

The method must be the same for the calibration batch and the current batch.

-Or-

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 from the shortcut menu.



 Table 55.
 Instrument states

| 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. | |
| Turn Device Off | Keeps the system in the Off state when the current run finishes. The Off state indicates that all power to the instrument, which the TraceFinder application can control, is turned off. This includes power to all heaters and subassemblies, but in some cases not all subassemblies. | |
| | 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. | |
| 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

From the Finish view, click Save.

The TraceFinder application saves your batch as a to-be-run (.tbr) file, closes the Acquisition mode, and returns you to the application dashboard.

To start an acquisition

1. Click Submit.

The Submit Options dialog box opens. For a detailed description of the parameters, see "Submit Options dialog box" on page 241.

2. (Optional) Select the **Create Reports** check box.

Note By default, the application acquires and processes data when you submit the batch.

- 3. Select the **Use** check box for the device that you want to use for this acquisition.
- 4. (Optional) Select the **Start Device** check box to indicate the device that will initiate communication with the other instruments.

This is usually the autosampler.

5. (Optional) 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.

6. (Optional with multiplexing enabled) 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 enabled) 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.
- 8. (Optional) Click Cancel to exit the Acquisition mode without performing any tasks.
- 9. To start the selected processes, click **OK**.

The selected processes begin, and the TraceFinder application returns you to the dashboard and shows the real-time display at the bottom of the dashboard.

The real-time display is visible from the dashboard and all modes. You can begin another batch in the Acquisition mode while you watch the real-time display of the currently acquiring batch.

Figure 67. Submit Options dialog box

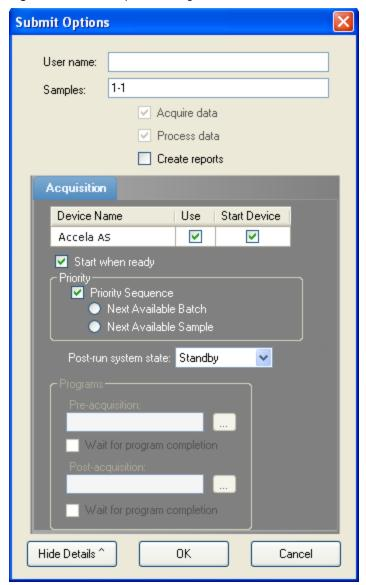


Table 56. Submit Options dialog box parameters (Sheet 1 of 2)

| Parameter | Description |
|----------------|---|
| User Name | Name of the current user. |
| Acquire Data | (Default) Submits the current batch to acquisition. |
| Process Data | (Default) Processes the data for the current batch. |
| Create Reports | Creates reports for the current batch. |

5 Using the Acquisition Mode Working with Batches

Table 56. Submit Options dialog box parameters (Sheet 2 of 2)

| Parameter | Description | |
|-----------------------|--|--|
| | Description | |
| Acquisition pane | | |
| Device Name | Lists all configured instruments. | |
| | If the instrument you want to use is not configured, close the TraceFinder application, configure the instrument, and then reopen the TraceFinder application. You cannot configure an instrument while the TraceFinder application is running. | |
| Use | Specifies the instruments used for this acquisition. | |
| Start Device | Specifies the instrument that initiates the communication with the other instruments. This is usually the autosampler. | |
| 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. | |
| Priority Sequence | With multiplexing enabled, places the batch immediately after the currently acquiring batch. | |
| | Without multiplexing enabled, specifies 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. | |
| Post-run system state | Specifies the system state after it acquires the last batch. On (default), Standby, or Off. | |
| 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

- The TraceFinder application writes saved batches to the subproject folder with the file extension .tbr (to be run):
 - ...\TraceFinder\1.1\Projects\projectname\subprojectname
- For each acquired sample, the application writes an RSX file to the batch Data folder:
 - $... \ \ | project name \ \ \ | batch name \ \ \ | Data$
- The application saves method information to the batch Methods folder:
 - $... \ \ | batchname \ \ | batchname \ \ | Methods \ | methodname$
- The application writes the reports to the batch Reports folder:
 - $... \ \ | project name \ \ | batch name \ \ | Reports$

5 Using the Acquisition Mode

Working with Batches

Figure 68. Finish view



Table 57. Finish view parameters (Sheet 1 of 2)

| Parameter | Description |
|---------------------------|--|
| System Status | The System Status pane displays the following: Devices used for the acquisition Project, subproject, and name of the batch Number of samples in the batch Number of standard and custom reports to be printed and saved as PDF, XML, or XLSM files Local method and instrument method used for the batch Number of compounds in the method |
| System Startup Method | The instrument method that runs before the batch. No autosampler injection takes place. This feature is not available for all instruments. |
| System Shutdown Method | The instrument method that runs after the batch. No autosampler injection takes place. This feature is not available for all instruments. |

Table 57. Finish view parameters (Sheet 2 of 2)

| Parameter | Description |
|-------------|--|
| Calibration | 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 to-be-run (.tbr) file. |
| Submit | Opens the Submit Options dialog box where you can optionally choose to generate reports. |

Real-time Display

You can access the real-time display from the dashboard and from any mode in the TraceFinder application.

❖ To access the real-time display from the dashboard

Click Real Time Status.



The real-time status displays at the bottom of the dashboard.

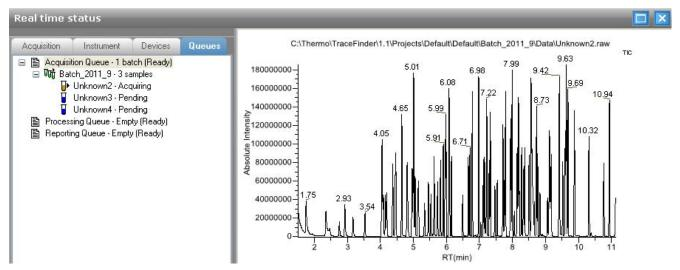
To access the real-time display from all modes

Click Real Time Status.



The real-time status displays at the bottom of the current view.

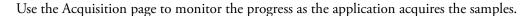
Figure 69. Real Time Status display

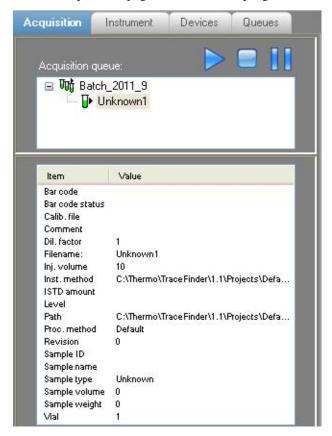


The real-time status display has four pages of information and a real-time trace pane:

- Acquisition Page
- Instrument Page
- Devices Page
- Queues Page
- Real-time Trace Display

Acquisition Page





❖ To pause or stop the batches in the queue

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

5 Using the Acquisition Mode

Real-time Display

Instrument Page

Use the Instrument page to monitor the currently acquiring sample.



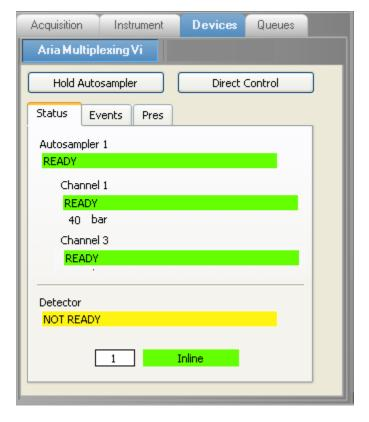
Devices Page

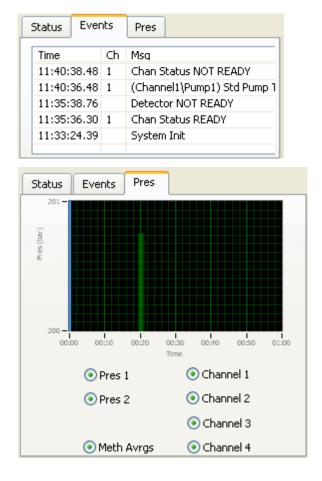
Use the Devices page to monitor the status of the instrument. The feedback you see on the Devices page depends on the instrument you are using. The following examples show an Accela autosampler and an Aria multiplexing device.

Accela Autosampler Feedback



Aria Multiplexing Feedback





5 Using the Acquisition Mode

Real-time Display

Use the following procedures:

- To pause the autosampler
- To access the Aria multiplexing controls
- To control the channels
- To view the pressure trace

To pause the autosampler

1. Click Hold Autosampler.

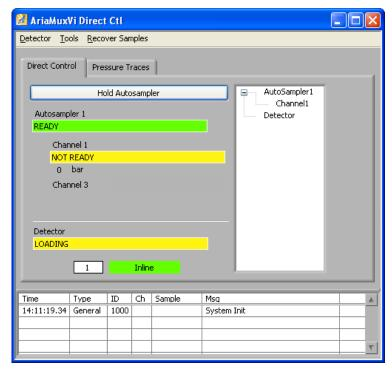
The autosampler finishes the current autosampler step and then pauses. The LC pumps and autosampler continue.

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

❖ To access the Aria multiplexing controls

Click Direct Control.

The AriaMuxVi Direct Ctl dialog box opens.



For a detailed description of the features in this dialog box, refer to the *Transcend Systems with Xcalibur Software User Guide*.

❖ To control the channels

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

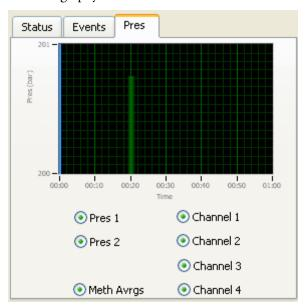


| Parameter | Description | |
|------------------|--|--|
| On | Turns on a stopped pump and continues acquiring the sample list assigned to that channel. | |
| Off | After the current sample completes, the application stops acquiring and the pump shuts down. | |
| Standby | After the current sample completes, the application stops acquiring. The pump continues to run. | |
| Disable / Enable | 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. Enable allows the channel to receive samples. | |
| | 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 . | |

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.

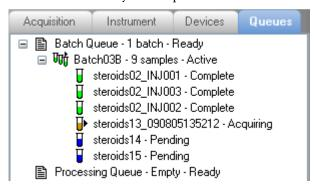


- To view the pressure for a specific pump, select the Pres 1 or Pres 2 option.
 By default, the pressure for all pumps are displayed.
- To view the pressure for a specific channel, select the corresponding channel number.By default, the pressure for all channels is displayed.

Queues Page

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

- Use the Queue-level Commands to pause or remove batches in any of the queues.
- Use the Batch-level Commands to pause or remove entire batches or samples within batches from any of the queues.



Queue-level Commands

Use the queue-level commands to pause or remove batches in any of the queues on the Queues page.

Use the following 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

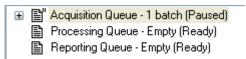
❖ To pause all batches in a queue

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

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

2. Right-click and choose **Pause Queue** from the shortcut menu.

After the current sample completes, 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** from the shortcut menu.

❖ To remove a single batch from a queue

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

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 completes, 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, Processing, or Reporting).
- 2. Right-click and choose **Stop All Batches** from the shortcut menu.

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, Processing, or Reporting).
- 2. Right-click and choose **Remove Pending Batches** from the shortcut menu.

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.

Figure 70. Queue-level shortcut menu

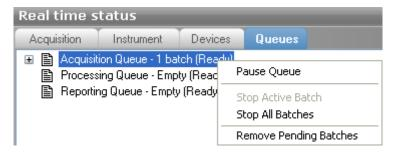


Table 58. Queue-level shortcut menu commands (Sheet 1 of 2)

| Parameter | Description |
|-------------------|--|
| Pause Queue | After the current sample completes, the application pauses the specified queue. Only the selected queue is affected. |
| Stop Active Batch | Removes all pending samples from the specified queue. The active sample is not affected. |

Table 58. Queue-level shortcut menu commands (Sheet 2 of 2)

| Parameter | Description |
|---------------------------|--|
| Stop All Batches | Removes all pending samples and batches from the specified queue. The active sample is not affected. |
| 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.

Use the following procedures:

- To stop a batch
- To remove a pending batch
- To remove pending samples from a batch
- To remove a single pending sample from a batch

To stop a batch

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

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

2. Right-click and choose **Stop Batch** from the shortcut menu.

The application confirms that you want to remove the selected batch from the queue. After the current sample completes, 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, 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 from the shortcut menu.

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

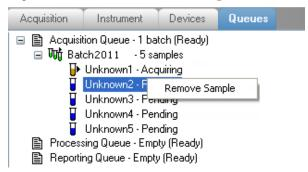
To remove pending samples from a batch

- 1. Select a batch in any of the queues (Acquisition, Processing, or Reporting).
 - The batch must have at least one pending sample.
- 2. Right-click and choose **Remove Pending Samples** from the shortcut menu.

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 remove a single pending sample from a batch

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



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

Figure 71. Batch-level shortcut menu

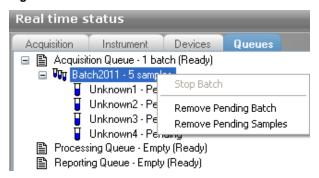
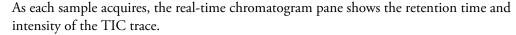
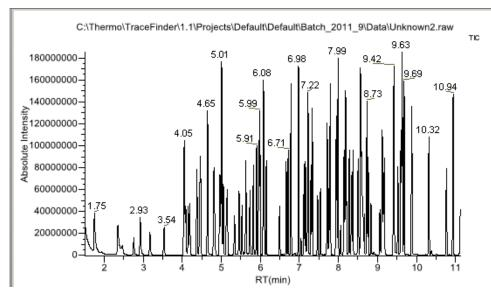


Table 59. Batch-level shortcut menu commands

| Parameter | Description |
|---------------------------|--|
| Stop Batch | After the current sample completes, 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. |

Real-time Trace Display



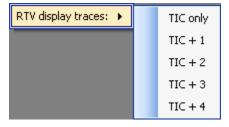


By default, the real-time display shows only the TIC trace as each sample acquires. To observe specific traces, such as the internal standard, use the RTV Display Traces function to display multiple traces.

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 Viewer" on page 156.

❖ To display multiple traces

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



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

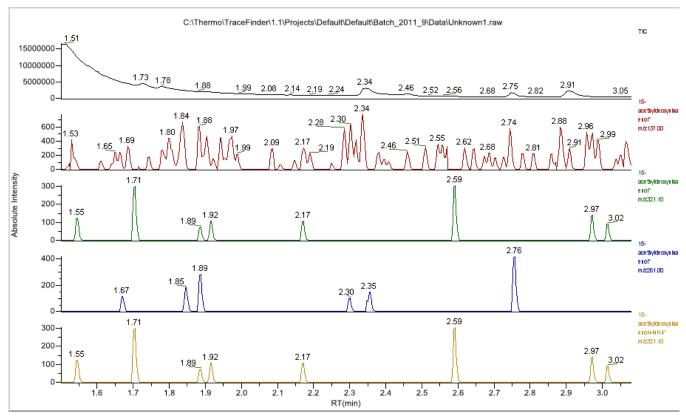
The TIC is always displayed at the top. When there are more traces than can fit in the pane, you can scroll through the traces.

For each trace, the application displays the mass or precursor mass.

5 Using the Acquisition Mode

Real-time Display

Figure 72. Real-time display with multiple traces



Sample Types

The TraceFinder application uses the following sample types in all sample definitions and reports. To view example standard reports specific to a sample type, see Appendix A, "Reports."

Table 60. Sample type definitions

| Sample type | Definition | |
|--------------|---|--|
| Matrix Blank | 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. | |
| Cal Std | (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 software can generate a calibration curve for each compound. | |
| Chk Std | (Check standard) 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 Chk Std sample, it compares the measured quantity with the expected value and an acceptability range. The quantitative analysis of a Chk Std sample is classified as <i>passed</i> if the difference between the observed and expected quantities is within the user-defined tolerance. A Chk Std 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. | |
| Unknown | Used for quantitative analysis of samples. | |
| Unknown/TIC | Used for quantitative and qualitative analysis of samples. | |
| LCS | Lab control sample. (Environmental and Food Safety only) | |
| LCSD | Lab control sample duplicate. (Environmental and Food Safety only) | |
| MDL | Method detection limits sample. (Environmental and Food Safety only) | |
| Method Val | Method validation sample. (Environmental and Food Safety only) | |
| MS | Matrix spike sample. (Environmental and Food Safety only) | |
| MSD | Matrix spike duplicate sample. (Environmental and Food Safety only) | |

Using the Data Review Mode

This chapter includes instructions for using the features of the Data Review mode.

Contents

- Working in the Batch View
- Working in Data Review View
- Working in Report View
- Working in the Local Method View

In the Data Review mode, you can review batches, batch data, reports, and local methods.

❖ To access the Data Review mode

Click **Data Review** from the dashboard or the navigation pane.



The Data Review navigation pane opens.



6 Using the Data Review Mode

Working in the Batch View

Working in the Batch View

In the Batch View, you can manually create and edit a new 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.

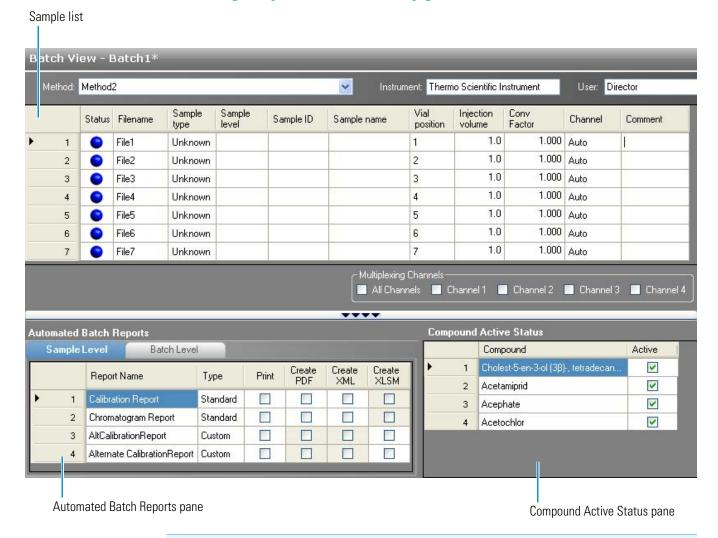
This section contains information about the following topics:

- Batch View Panes
- Creating a New Batch
- Editing a Batch
- Editing Report Output Formats
- Submitting a Batch

Batch View Panes

The Batch View is divided into three panes:

- Use the sample list to create a batch. See "Creating a New Batch" on page 264.
- Use the Automated Batch Reports pane to select the type of output formats you want to generate for the reports. See "Editing Report Output Formats" on page 274.
- Use the Compound Active Status pane to make specific compounds active or inactive. See "Setting Compound Active Status" on page 276.



Tip To resize the panes, drag the separators that divide the panes.

Creating a New Batch

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

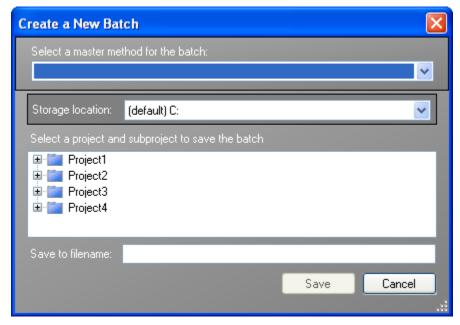
Use the following procedures:

- 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 customize the column display
- To temporarily customize the column display

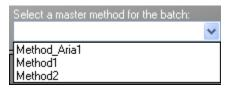
❖ To create a new batch

1. Click **New Batch** in the Batch View task pane or choose **File > New > Batch**.

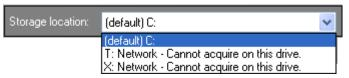
The Create a New Batch dialog box opens.



2. Select a master method from the Method list.



3. Select a drive from the Storage Location list.



The project list displays all projects, subprojects, and batches on the selected drive.

Tip The application does not display drives that do not have a project and subproject.

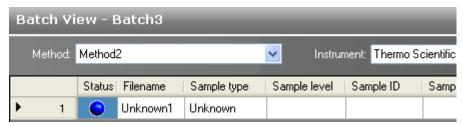
You cannot use network drives to acquire data. For more information about network drives, see "Working with Drives" on page 40.

4. Select a project and a subproject and enter a name for your new batch.

Tip To enable the Save button, you must select a subproject and enter a unique batch name. If the Save button is not enabled, either you have entered a name that is already used or you have not selected a subproject.

5. Click Save.

A new, unnamed batch opens with one Unknown sample.



❖ To add samples to the list

- 1. To add a single sample row, right-click the sample list and choose **Add Sample** from the shortcut menu.
- 2. To add multiple sample rows, select the number of rows and then click the **Add Sample** button, 3 .

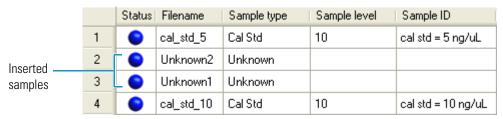
The application adds the specified number of new, empty samples to the end of the sample list.

❖ To insert samples into the list

Select the sample above which you want to insert unknown samples, and do one of the following:

- To insert a single sample row, right-click and choose **Insert Sample** from the shortcut menu.
- To insert multiple sample rows, select the number of rows and then click the **Insert Sample** button 3 .

The TraceFinder application inserts a new, empty sample or samples above the selected sample.

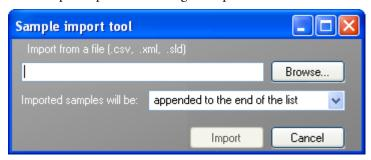


❖ To import samples into the list

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



The Sample Import Tool dialog box opens.



From this dialog box, you can import samples from a .csv, .xml, or .sld file.

- Click Browse and select a .csv, .xml, or .sld file that contains the samples you want 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 samples list.

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

| Xcalibur column | TraceFinder column |
|-----------------|--------------------|
| Level | Sample Level |
| Position | Vial Position |
| Inj Vol | Injection Volume |
| Dil Factor | Conv Factor |

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

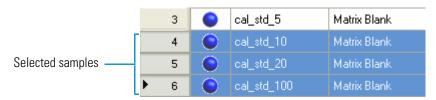
| Xcalibur sample type | TraceFinder sample type |
|----------------------|-------------------------|
| Blank | Matrix Blank |
| QC | Chk Std |
| Std Bracket | Cal Std |

(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 you want to remove.

Make sure the first column indicates that the samples are selected.



2. Right-click and choose **Remove Selected Samples** from the shortcut menu.

❖ To copy a sample

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

 The TraceFinder application inserts the copy above the selected sample.

To reinject a sample

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

The TraceFinder 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 TraceFinder application copies all parameter values from the original sample.

❖ To edit sample values

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

Double-click the Filename column and type a file name.

-Or-

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

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

| Available Environmental and Food Safety sample types | | | | | |
|--|-------------|-------------|--|--|--|
| Matrix Blank | Solvent | Unknown/TIC | | | |
| Cal Std | Chk Std | Unknown | | | |
| LCS | MDL | MS | | | |
| LCSD | Method Val | MSD | | | |
| Available General Quantitation sample types | | | | | |
| Matrix Blank | Unknown/TIC | Unknown | | | |
| Cal Std | Chk Std | Solvent | | | |

3. For each Cal Std or Chk Std sample, click the Sample Level cell and select a level from the list.

The sample levels are defined in the master method. If there is nothing to select in the Sample Level list, do the following:

- a. Return to the Method Development mode.
- b. Open the method.
- c. Click the Calibration Levels tab.
- d. Add the levels.
- e. Save the method.
- f. Return to the Data Review mode, and click **Update**.

For detailed instructions, see Chapter 4, "Using the Method Development Mode."

4. Enter or edit the values for the remaining columns.

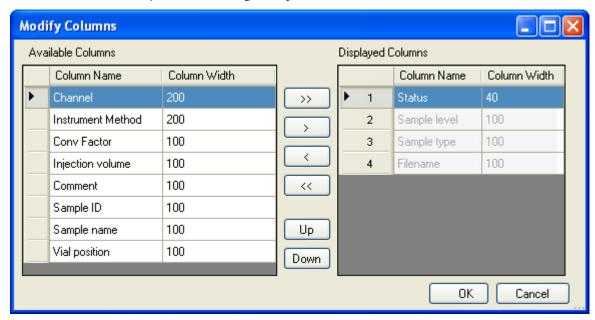
When you use the scroll bar at the bottom of the samples list, the Status, Flags, Filename, and Sample Type columns remain fixed and only the other columns scroll right and left.

For instructions to automatically copy or fill values in these columns, see Appendix B, "Using Copy Down and Fill Down."

To customize the column display

1. Right-click the Batch View sample list and choose **Modify Columns** from the shortcut menu.

The Modify Columns dialog box opens.



| Parameter | Description | |
|---------------|--|--|
| >> | Moves all columns to the Displayed Columns pane. | |
| \rightarrow | Moves the selected column to the Displayed Columns pane. | |
| < | Moves the selected column to the Available Columns pane. You cannot move the Status, Sample Level, Sample Type, or Filename columns. | |
| << | Moves all columns except Status, Sample Level, Sample Type, and Filename to the Available Columns pane. | |
| Up | Moves the selected column name in the Displayed Columns pane one row up in the column order. You cannot move the Status, Sample Level, Sample Type, or Filename columns. | |
| Down | Moves the selected column name in the Displayed Columns pane one row down in the column order. You cannot move the Status, Sample Level, Sample Type, or Filename columns. | |

6 Using the Data Review Mode

Working in the Batch View

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

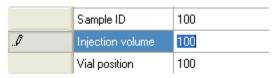
All the columns you select are displayed after the Status, Sample Level, Sample Type, and Filename columns.

- 3. To arrange the order of the columns, do the following:
 - a. In the Displayed Columns pane, select a column name.
 - b. Use the **Up** and **Down** buttons 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; the last column in the list represents the rightmost column in the Batch View sample list.

Note You cannot move the Status, Sample Level, Sample Type, or Filename columns.

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



- 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.

You can use the drag-and-drop method to temporarily reorder the columns in your sample list, but you cannot save this order. When you restart the TraceFinder application, the Batch View displays your columns in the order specified in the Modify Columns dialog box.

❖ To temporarily customize the column display

- 1. To resize a column, drag the header separator on the right side of the column.
- 2. To move a column, drag the column header.

You can swap the Status, Sample Level, Sample Type, and Filename columns with each other, but they must always be the first four columns in the sample list.

You cannot save these changes. When you restart the TraceFinder application, the Batch View displays your columns in the order specified in the Modify Columns dialog box.

Editing a Batch

In the Batch View, you can open a saved batch and edit the samples 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 to-be-run (.tbr) batch.

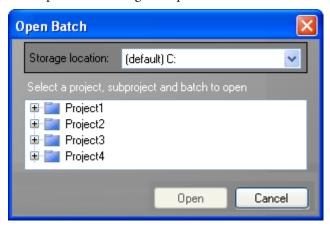
Use the following procedures:

- 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 submit all samples in the batch
- To submit selected samples

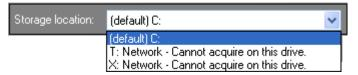
To open a saved batch

1. From the Batch View task pane, click Open Batch.

The Open Batch dialog box opens.



2. Select a drive from the Storage Location list.



The project list displays all projects, subprojects, and batches on the selected drive.

- 3. Select a project, subproject, and batch.
- 4. Click Open.

The selected batch opens in the Batch View.

❖ To open a recent batch

Click a batch name in the Recent Files list.

The selected batch opens in the Batch View.

Tip To view the drive, project, and subproject for a recent batch, hold your cursor over the batch name.

Batch View

New batch

Open batch

Recent files

Batch3

Drive: C:
Project: Project1
Subproject: SubProject1
Batch: Batch1

❖ To edit samples in a batch

Use the commands described in "Working in the Batch View" on page 262.

You can edit samples, add new samples, reinject acquired samples, or delete samples.

To reinject a sample from a previously acquired batch

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

The TraceFinder 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 TraceFinder 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_std_50_INJ001 | Cal Std | 10 |
|---|--------------------|---------|----|
| | cal_std_50_INJ002 | Cal Std | 10 |
| • | cal_std_50 | Cal Std | 10 |
| | cal_std_100_INJ001 | Cal Std | 10 |
| • | cal_std_100 | Cal Std | 10 |

When you submit all samples in this batch, all samples (including previously acquired samples) are acquired.

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 .tbr batch. You can open this batch from the Ready to Acquire page in the Acquisition mode and submit the batch. All submitted samples—both the reinjection samples and the previously acquired samples—are acquired. The application appends a timestamp to the acquired raw data files to differentiate each acquisition.

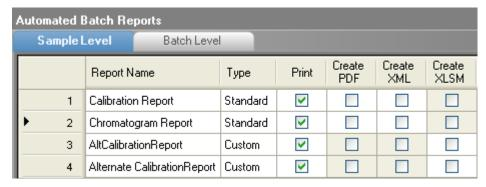
Editing Report Output Formats

In the Automated Batch Reports pane, you can view the reports that are selected for this batch and modify which output formats are generated for each report.

❖ To edit the sample-level output formats

1. Click the **Sample Level** tab.

The application displays reports and the output formats as they were specified in the method.



For detailed instructions for specifying which reports and output formats are generated, see "Specifying the Reports Configuration" on page 58.

- 2. Select or clear any of the check boxes for your reports.
- 3. To duplicate an output format for all reports for this sample, right-click the cell and choose **Copy Down** from the shortcut menu.

All check boxes in the column below the selected cell duplicate the selected or cleared state in the selected cell. You can duplicate the output type only for reports that have this output format available.

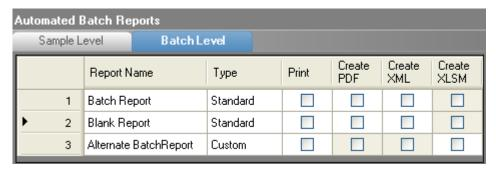
4. To duplicate the output format for all samples in the batch, right-click the cell and choose **Apply Selection to All Samples** from the shortcut menu.

You can duplicate the output type only for reports that have this output format available.

❖ To edit the batch-level output formats

1. Click the Batch Level tab.

The application displays the reports and the output formats as they were specified in the method.



For detailed instructions for specifying which reports and output formats are generated and which reports are batch-level, see "Specifying the Reports Configuration" on page 58.

- 2. Select or clear any of the check boxes for your reports.
- 3. To duplicate the output format for all reports, right-click the cell and choose **Copy Down** from the shortcut menu.

All check boxes in the column below the selected cell duplicate the selected or cleared state in the selected cell. You can duplicate the output type only for reports that have this output format available.

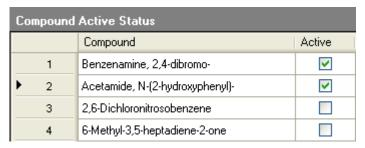
Setting Compound Active Status

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

❖ To set a compound as active or inactive

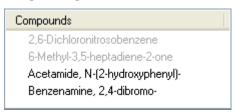
1. In the samples list, select a sample.

All compounds in the selected sample are listed in the Compound Active Status pane. The default active/inactive status is determined by the settings in the local method. See "Identification" on page 119.



2. Select or clear the Active check box for the compound.

When you specify a compound as inactive in this pane, it becomes inactive in the Compounds pane in the Data Review view. Inactive compounds are grayed out:



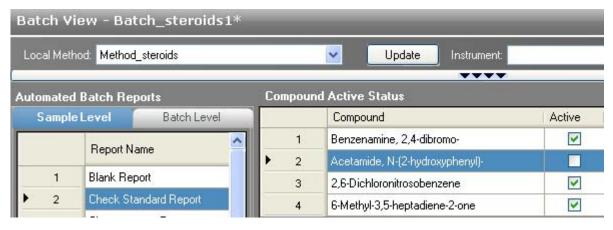
For instructions for changing the active/inactive status in the Data Review view, see "Inactive and Excluded Compounds" on page 290.

Compound Active/Inactive Status

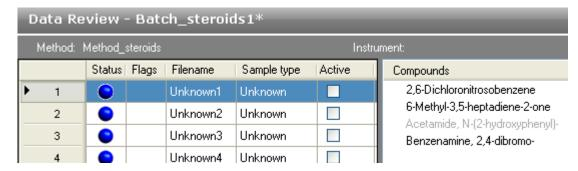
You can specify which compounds are active or inactive in the Local Method View, the Batch View, or the Data Review view.



For details about the setting the status on the Identification page, see "Identification" on page 119.



For details about setting the status in the Batch View, see "Setting Compound Active Status" on page 276.



For details about setting the status in the Data Review view, see "Inactive and Excluded Compounds" on page 290.

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" on page 280.

Use the following procedures:

- To submit all samples in the batch
- To submit selected samples

For a description of commands on the shortcut menu, see "Batch View shortcut menu commands" on page 282.

❖ To submit all samples in the batch

1. Click the **Submit Batch** button,

The Submit Options dialog box opens. See "Submit Options dialog box" on page 280.

- 2. Select the tasks you want to perform: acquire data, process data, or create reports.
- 3. (Optional) Click **Show Details** to display additional Acquisition parameters.
- 4. Select the **Use** check box for the device that you want to use for this acquisition.
- 5. (Optional) Select the **Start Device** check box to indicate the device that will initiate the communication with the other instruments.

This is usually the autosampler.

(Optional) Select the **Start When Ready** check box to have all instruments start together when they are all ready.

When this is cleared, individual instruments can start at different times and then must wait for the last instrument to be ready.

7. (Optional with multiplexing enabled) Select the **Priority Sequence** check box.

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

- 8. (Optional without multiplexing enabled) 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.
- 9. To start the selected processes, click **OK**.

To submit selected samples

- 1. Select the samples you want to submit.
- Click the Submit Selected Samples button, The Submit Options dialog box opens. See "Submit Options dialog box" on page 280.
- 3. Select the tasks you want to perform: acquire data, process data, or create reports.
- 4. Select the **Use** check box for the device that you want to use for this acquisition.
- 5. (Optional) Click **Show Details** to display additional Acquisition parameters.
- 6. (Optional) Select the **Start Device** check box to indicate the device that will initiate communication with the other instruments.
 - This is usually the autosampler.
- 7. (Optional) Select the **Start When Ready** check box to have all instruments start together when they are all ready.
 - When this is cleared, individual instruments can start at different times and then must wait for the last instrument to be ready.
- 8. (Optional with multiplexing enabled) Select the **Priority Sequence** check box.
 - The application acquires the priority batch on the next available channel or the assigned channel.
- 9. (Optional without multiplexing enabled) 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.
- 10. To start the selected processes, click **OK**.

Submit Options User name: 1-1 Samples: Acquire data ✓ Process data Create reports Acquisition Device Name Use Start Device V V Accela AS Start when ready Priority-Priority Sequence Next Available Batch Next Available Sample Post-run system state: Standby Hide Details ^ OΚ Cancel

Figure 73. Submit Options dialog box

Table 61. Submit Options dialog box parameters (Sheet 1 of 2)

| Parameter | Description |
|----------------|---|
| User Name | Name of the current user. |
| Samples | Reports the batch number and number of samples. |
| Acquire Data | (Default) Submits the current batch to acquisition. |
| Process Data | (Default) Processes the data for the current batch. |
| Create Reports | Creates reports for the current batch. |

Table 61. Submit Options dialog box parameters (Sheet 2 of 2)

| Parameter | Description |
|-----------------------|--|
| Acquisition pane | |
| Device Name | Lists all configured instruments. |
| | If the instrument 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. |
| Use | Specifies the instruments used for this acquisition. |
| Start Device | Specifies the instrument that will initiate communication with the other instruments. This is usually the autosampler. |
| 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. |
| Priority Sequence | With multiplexing enabled, places the batch immediately after the currently acquiring batch. |
| | Without multiplexing enabled, specifies 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. |
| 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. |

6 Using the Data Review Mode

Working in the Batch View

Figure 74. Batch View

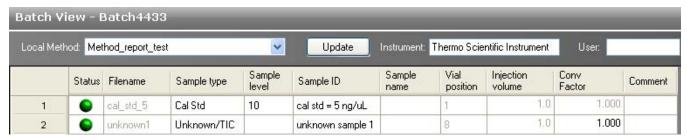


Table 62. Batch View shortcut menu commands (Sheet 1 of 2)

| 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" on page 266. |
| Browse in raw file | Opens a dialog box where you can select a raw data file to use for the sample row. |
| Map raw files to samples | Opens a dialog box where you can select raw data files to use for the 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. See "Column Display" on page 286. |
| Сору | Copies the data in the selected rows or columns to the Clipboard. Use this command to copy sample information to another application, such as an Excel spreadsheet. You cannot paste this data back into the Batch View sample list. |

Table 62. Batch View shortcut menu commands (Sheet 2 of 2)

| Command | Description | | |
|--------------------|---|--|--|
| 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 another application, such as an Excel spreadsheet. You cannot paste this data back into the Batch View sample list. For example: | | |
| | Sample type | | |
| | Matrix Blank | | |
| | Cal Std | | |
| | Chk Std Sample type | | |
| | Unknown/TIC Unknown/TIC | | |
| | Unknown/TIC | | |
| | Copy With Headers from TraceFinder Paste into Excel spreadsheet | | |
| Paste | Pastes a single column of copied data from another application, such as an Excel spreadsheet, 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. | | |

Working in Data Review View

In the Data Review view, you can view the data generated by the master method. Use the Data Review view to verify the data for a sample-specific compound before you generate reports. You can use the functions in the Data Review view to investigate and edit the quantification values in a batch.

❖ To open the Data Review view

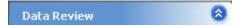
1. Do one of the following (to open Data Review mode):

From the dashboard, click Data Review.

-Or-

Click **Data Review** in the navigation pane of the current mode.

2. In the Data Review navigation pane, click **Data Review** (to open the Data Review view).



The Data Review view for the currently selected batch opens.

The Data Review view uses a samples list and one of two modes: Quan Mode or Qual Mode. The Qual Mode is available only for Unknown/TIC sample types. When you view the data for an Unknown/TIC sample type, you can switch between Qual Mode and Quan Mode.

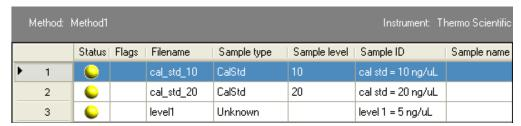
This section includes the following topics:

- Samples List
- Quan Mode
- Qual Mode

Samples List

Use the samples list to select a particular sample. For detailed descriptions of the columns in a samples list, see "Samples list" on page 292.

Status indicators for each sample indicate if the sample is unacquired, acquired, currently acquiring, or processed.



The samples list includes the following features:

- Column Display
- Status Indicators
- Sample Flags
- Compound Flags
- Inactive and Excluded Compounds

The samples list is the same in both Quan Mode and Qual Mode and displays all the quantitative data for the samples of a batch.

• In Quan Mode, the samples list works with the Compounds pane to select a unique sample and compound combination, which then has its textual and graphical values displayed in the Quan Mode pane. The list of compounds that are available for a specific method is displayed in the Compounds pane.

From the samples list, you can make a compound active or inactive. Switching a compound to inactive status does not remove its data and calculated values from the result set. Instead, the TraceFinder application masks the appearance of that compound for that particular sample and grays the compound in the Compounds list. For a calibration standard, 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. It is no longer part of the result set.

In Qual Mode, the samples list works with the peaks pane to select a unique sample and
peak combination, which then has its textual and graphical values displayed in the Qual
Mode pane.

6 Using the Data Review Mode

Working in Data Review View

Column Display

The samples list can contain many columns of information. You can scroll to see all the columns of information, and you can customize which columns you want to display and their display order.

Use the following procedures:

- To scroll the samples list
- Status Indicators
- To customize the column display

❖ To scroll the samples list

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

When you use the horizontal scroll bar at the bottom of the samples list, the Status and Filename columns stay fixed while the other columns scroll right and left.

To temporarily customize the column display

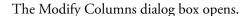
- 1. To resize a column, drag the header separator on the right side of the column.
- 2. To move a column, drag the column header.

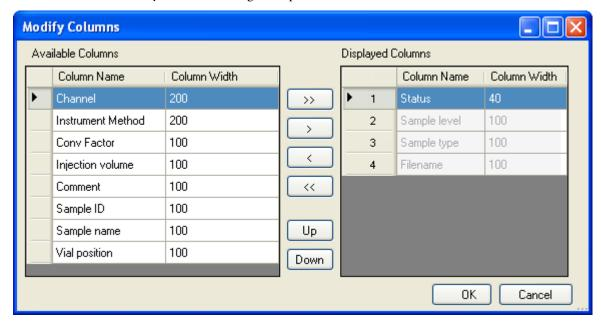
The Status, Sample Level, Sample Type, and Filename columns can be swapped with each other, but they must always be the first four columns in the sample list.

You cannot save these changes. When you restart the TraceFinder application, the Data Review view displays your columns in the order specified in the Modify Columns dialog box.

❖ To customize the column display

1. Right-click the Data Review sample list and choose **Modify Columns** from the shortcut menu.





| Parameter | Description |
|-----------|--|
| >> | Moves all columns to the Displayed Columns pane. |
| > | Moves the selected column to the Displayed Columns pane. |
| < | Moves the selected column to the Available Columns pane. You cannot move the Status, Sample Level, Sample Type, or Filename columns. |
| << | Moves all columns except Status, Sample Level, Sample Type, and Filename to the Available Columns pane. |
| Up | Moves the selected column name in the Displayed Columns pane one row up in the column order. You cannot move the Status, Sample Level, Sample Type, or Filename columns. |
| Down | Moves the selected column name in the Displayed Columns pane one row down in the column order. You cannot move the Status, Sample Level, Sample Type, or Filename columns. |

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

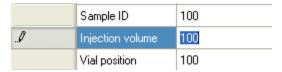
All the columns you select are displayed after the Status, Sample Level, Sample Type, and Filename columns.

- 3. To arrange the order of the columns, do the following:
 - a. In the Displayed Columns pane, select a column name.
 - b. Use the **Up** and **Down** buttons 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; the last column in the list represents the rightmost column in the Batch View sample list.

Note You cannot move the Status, Sample Level, Sample Type, or Filename columns.

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



- b. Type a new value for the width.
- 5. When you have completed your changes, click **OK**.

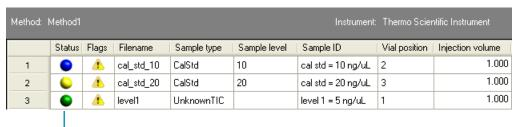
The columns in the sample list immediately reflect your changes. The application uses these settings for all sample lists in the Data Review view.

You can use the drag-and-drop method to temporarily reorder the columns in your sample list, but you cannot save this order. When you restart the TraceFinder application, the Data Review view displays your columns in the order specified in the Modify Columns dialog box.

Status Indicators

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

- · Orange for samples that are currently acquiring
- Blue for samples that have not been acquired
- Yellow for samples that have been acquired and not processed
- Green for samples that have been acquired and processed

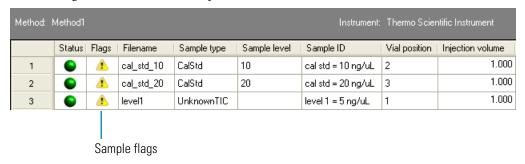


Status indicators

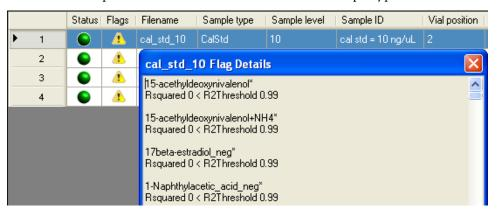
Sample Flags

The Flags column in the samples list displays a caution flag if the sample is not in compliance with the method criteria.

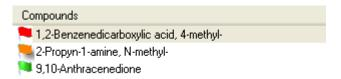
Sample caution flags remain static when you switch between compounds for chromatogram review until a change is completed, for example, when a compound is manually integrated and no longer falls outside the accepted criteria.



To display a summary of problems found in the sample, click the caution flag. The summary does not list compounds that are not found in Unknown sample types.



Compound Flags



Flags are displayed in these situations:

- When a compound has violated (or is activated by) any of the values set in the method (See "Editing the Flags Page" on page 159.)
- For compounds that are not found
- For compounds that are not found in Cal Std or Chk Std sample types
- For compounds that are outside the specified ion ratio range

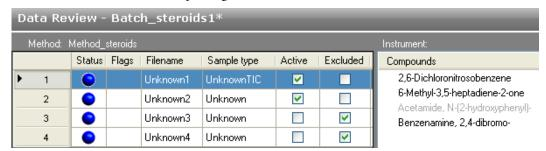
Working in Data Review View

The compounds list is sorted first by flag indicators and then by compound names. Compound flags indicate the following:

- Red flags for compounds that have ion ratio failures or method validation failures
- Orange flags for compounds that are below the LOQ, below the LOD, or between the LOD and LOQ values specified in the method
- Green flags for compounds that are over the LOR amount specified in the method
- No flag for compounds that have no errors or where no report options are selected

Inactive and Excluded Compounds

Use the Active and Excluded columns to control which compounds are used for calculating the calibration curve and for reporting.



Use the following procedures:

- To make a sample active or inactive
- To exclude a calibration point

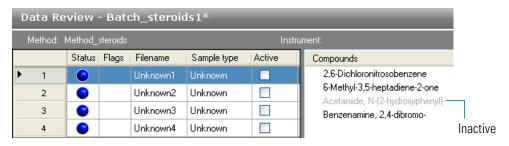
To make a sample active or inactive

1. Select the sample in the samples list.

All compounds in the selected sample are displayed 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.
- 3. In the samples list, 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 samples list, select the **Excluded** check box for the sample.

Use the horizontal scroll bar at the bottom of the table to scroll to the Excluded column.

When a value is no longer used for calibration, it is displayed as an empty box in the graphical view of the calibration curve.

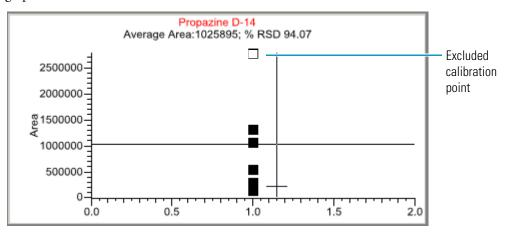


Figure 75. Samples list

| Method | Method: Batch1680B_Steroid Method | | | | | Instru | ment: Therm | o Scientifi | c Instrument | |
|--------|-----------------------------------|----------|------------|-------------|-----------------|--------------|----------------|-------------|------------------|---------------------|
| | Status | Flags | Filename | Sample type | Sample level | Sample ID | Sample name | Comment | Vial position | Injection volume |
| 1 | • | <u> </u> | steroids02 | Unknown | | Sample02 | | | 1 | 20.0 |
| 2 | • | <u> </u> | steroids03 | Unknown | | Sample03 | | | 2 | 20.0 |
| 3 | • | | steroids04 | Unknown | | Sample04 | | | 3 | 20.0 |
| 4 | • | | steroids05 | Unknown | | Sample05 | | | 4 | 20.0 |

| Integration mode | Height | *Area | Actual RT | Expected RT | Calc Amt | Theo Amt | Resp ratio | IS Amt | IS Resp |
|---------------------|--------|---------|--------------|----------------|----------|----------|---------------|--------|------------|
| Method | 312402 | 2958671 | 3.16 | 3.19 | N/A | | 4.372 | 0.600 | 676792 |
| Method | 152082 | 1590374 | 3.17 | 3.19 | N/A | | 2.361 | 0.600 | 673528 |
| Method | 71110 | 766235 | 3.15 | 3.19 | N/A | | 0.992 | 0.600 | 772347 |
| Method | 583384 | 6056721 | 3.11 | 3.19 | N/A | | 9.472 | 0.600 | 639426 |

| Active | Excluded | % Diff | % RSD | % CV | Channel |
|----------|----------|--------|-------|------|-----------|
| ~ | | N/A | N/A | N/A | Channel 1 |
| ✓ | | N/A | N/A | N/A | Channel 3 |
| ✓ | | N/A | N/A | N/A | Channel 1 |
| ✓ | | N/A | N/A | N/A | Channel 3 |

Cells in the samples list that should not have a value, such as the theoretical concentration for an unknown, are shaded and empty. Cells that should have a value, but none exists, report N/A (not available). Results for compounds that are not detected display N/F (not found).

Table 63. Samples list columns (Sheet 1 of 3)

| Parameter | Description |
|-------------|--|
| Status | Orange for samples that are currently acquiring Blue for samples that have not been acquired Yellow for samples that have been acquired and not processed Green for samples that have been acquired and processed |
| Flags | Displayed only when a compound within the sample has an error. |
| Filename | Name of the raw data file that contains the sample data. |
| Sample Type | Defines how the TraceFinder application processes the sample data. Each sample is classified as one of the following sample types: |
| | Environmental and Food Safety: Matrix Blank, Cal Std, Chk Std, LCS, LCSD, MDL, Method Val, MS, MSD, Unknown, Unknown/TIC, Solvent |
| | General Quantitation: Matrix Blank, Cal Std, Chk Std, Unknown, Unknown/TIC, Solvent |

Table 63. Samples list columns (Sheet 2 of 3)

| Parameter | Description |
|------------------|---|
| Sample 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. |
| Comment | A user-defined comment for the sample. |
| Vial Position | The tray vial number used for the autosampler acquisition. |
| Injection Volume | The injection volume in microliters of sample to be injected. |
| | 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 range displayed. The Injection Volume value set in the master method overwrites the value in the instrument method. |
| | Range: 0.1 through 5000 μL |
| Integration Mode | Indicates whether the peaks have been manually integrated or integrated from the original method. |
| Height | The distance from the peak maximum to the peak base, measured perpendicular to the ordinate. When the Resp Ratio is specified as Height, this column displays an asterisk (*Height). |
| Area | The area obtained by integrating peak intensities from the start to the end of the peak. When the Resp Ratio is specified as Area, this column displays an asterisk (*Area). |
| 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. |
| Expected RT | Expected retention time for the compound. |
| Calc Amt | The amount present in the sample, as determined using the calibration curve and the response ratio. |
| Theo Amt | Theoretical amount of the compound expected in the sample. |
| Resp 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. |
| IS Amt | Amount of internal standard. |
| 10 1 11111 | |

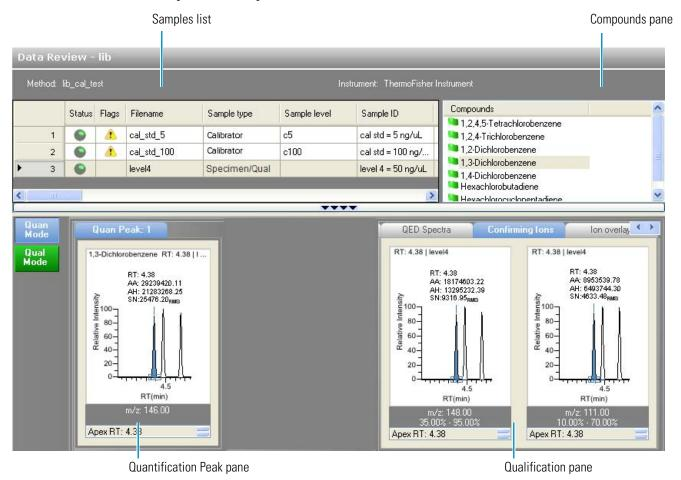
6 Using the Data Review Mode Working in Data Review View

Table 63. Samples list columns (Sheet 3 of 3)

| Parameter | Description |
|-----------|---|
| Active | Displays or hides a compound for a particular sample. 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. It is no longer part of the result set. |
| Excluded | Turns a compound on or off in the Calibration curve of the Qualification pane. |
| %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. |
| %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. |
| Channel | Specifies the channel on which the sample was run. If the sample is not acquired, the value is Pending. This column is displayed only when you enable multiplexing. |

Quan Mode

Use the Quan Mode and the associated Compounds pane to view quantitative information to complement the textual information for the selected sample. The Quan Mode displays quantitative peak and confirming ion information for selected compounds that are found in the processed samples.



In addition to the Samples list, the Quan Mode view uses the following panes:

- Compounds
- Quantification Peak
- Qualification

6 Using the Data Review Mode

Working in Data Review View

Compounds

The Compounds pane works with the samples list pane to display textual and graphical values for a unique file and compound combination.

Use the following procedures:

- To sort the compounds list
- To display peaks for a specific compound
- To display specific problems with a compound
- To display the internal standard for a compound

❖ To sort the compounds list

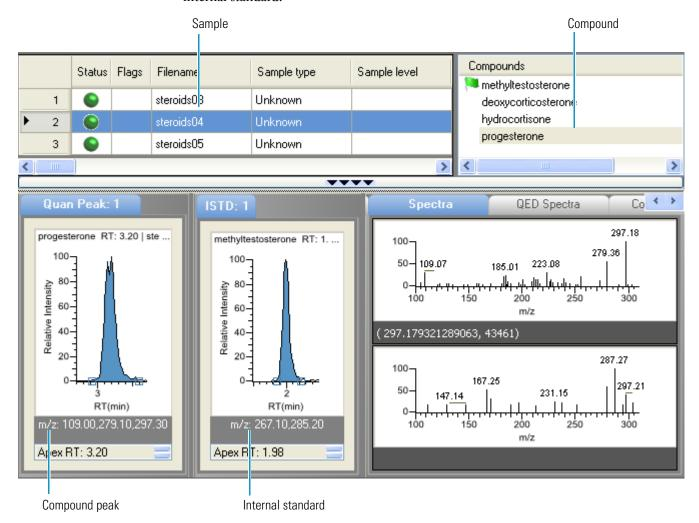
In the Compounds pane, right-click and choose one of the following sort styles from the shortcut menu:

| Command | Description |
|---------------------------------|---|
| Sort by flag and alphabetical | Sorts the compounds first by flag and then within each flag group, sorts the compounds alphabetically. |
| Sort by flag and retention time | Sorts the compounds first by flag and then within each flag group, sorts the compounds by retention time. |
| Sort by alphabetical | Sorts the compounds alphabetically $(1-n \text{ followed by } a-z)$. |
| Sort by retention time | Sorts the compounds from shorter retention time to longer retention time. |

To display peaks for a specific compound

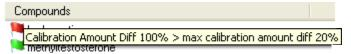
- 1. In the samples list, select the sample.
 - The Compounds pane lists all compounds specified in the method.
- 2. In the Compounds pane, select the compound in the sample.

The Quantification Peak pane displays the peaks for the selected compound and its internal standard.



To display specific problems with a compound

Hold the cursor over the flag to display the problems with the compound.



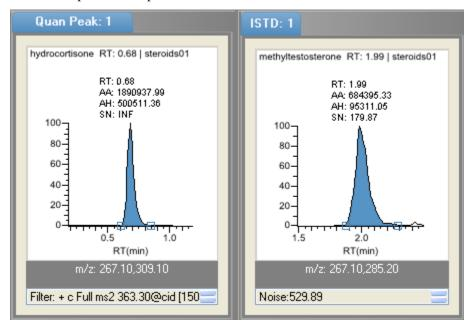
❖ To display the internal standard for a compound

1. Right-click the Quan Peak pane and choose **Show Internal Standard** from the shortcut menu.



Note By default, the ISTD pane is not displayed.

The Quantification Peak pane displays an ISTD pane with the internal standard for the selected sample and compound.



2. To hide the ISTD pane, right-click the Quan Peak or ISTD pane and choose **Hide Internal Standard** from the shortcut menu.

Quantification Peak

The Quantification Peak pane displays the compound selected in the samples list and Compound pane. 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 method values were originally calculated based on the processing method parameters. The manual values are a result of what has been edited.

When the sample contains an internal standard, the chromatogram shows both the analyte and the internal standard in side-by-side panes.

Use the following procedures:

- To zoom in on a peak
- To manually integrate a quantification or qualification ion
- To manually add a peak
- To remove a manually created peak
- To switch between method and manual integration modes
- To change the displayed information for detected peaks
- To modify the peak detection settings

For a description of commands on the shortcut menu, see "Quantification Peak Shortcut Menu" on page 303.

To zoom in on a peak

- In the chromatogram plot, drag your 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** from the shortcut menu.

To manually integrate a quantification or qualification ion

1. Place your cursor over one of the two peak delimiter tags in the Quantification 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.

Both the Quantification Peak pane and the Integration mode column in the Quantification Data pane reflect the change between method and manual modes. The generated reports for these data identify the manual modifications.

Note Because a Blank Report displays only the quan mass, when you manually integrate a confirming ion, the manual integration flag in the report is displayed on the quan mass.

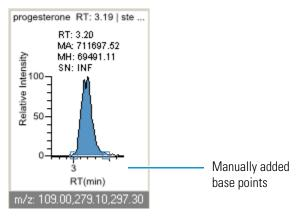
❖ To manually add a peak

1. Right-click anywhere in the Quantification Peak pane, and choose **Add Peak** from the shortcut menu.

If a peak is already detected, the Add Peak command is not enabled.

2. Click to indicate the left and right base points for the peak.

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 remove a manually created peak

Right-click the chromatogram plot, and choose **Remove Peak** from the shortcut menu.

The application removes the manually added peak.

To switch between method and manual integration modes

Right-click the chromatogram view and choose **Method Integration Settings** or **Manual Integration Settings** from the shortcut menu.

Initially, the method 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, the 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 these data identify the manual modifications.

Note Because a Blank Report displays only the quan mass, when you manually integrate a confirming ion, the manual integration flag in the report is displayed on the quan mass.

To change the displayed information for detected peaks

- 1. Right-click the quantification chromatogram plot and hold the cursor over Peak Labels.
- 2. Choose to display labels for the peak retention time, peak height, peak area, or signal to noise.

The label types in the list are selected for displayed labels and cleared for labels that are not displayed.

3. To remove a label, select the label type again and clear it.

Label settings are globally applied to quantification peaks, confirming peaks, and internal standard peaks.

Tip The labels do not always update on all peak displays. To update all labels, select a different compound, and then reselect the compound whose labels you changed.

To modify the peak detection settings

1. Right-click the chromatogram view and choose **Peak Detection Settings** from the shortcut menu.

The Peak Detection Settings dialog box opens. Use the Peak Detection Settings dialog box to adjust detection settings that were specified in the method.

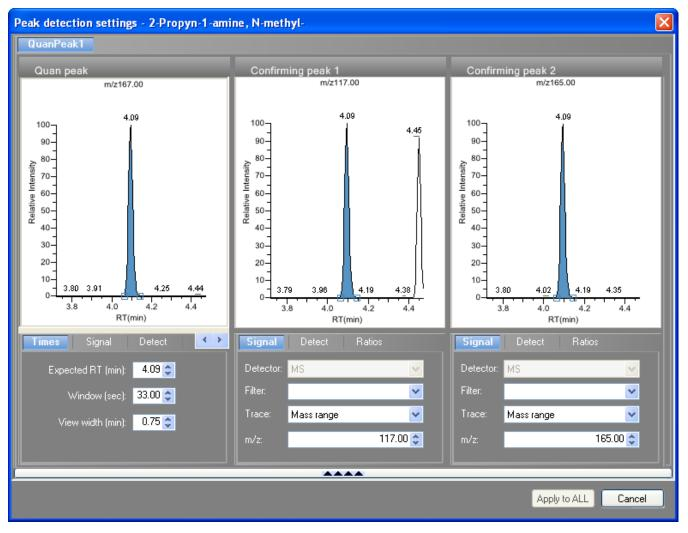


Figure 76. Peak Detection Settings dialog box

- Edit any of the detection settings.
 For detailed descriptions of all detection settings, see "Detection" on page 121.
- 3. To save your changes to this compound in all samples in this batch, click **Apply to All**.

Quantification Peak Shortcut Menu

The Quantification Peak shortcut menu includes the following commands:

Table 64. Quantification Peak shortcut menu commands

| Command | Description |
|--|---|
| Method Integration Settings | Displays method integration settings. |
| Manual Integration Settings | Displays manual integration settings. |
| Add Peak -or- Remove Peak -or- Cancel Add Peak | Adds a peak, removes a peak, or cancels an add peak operation in progress. |
| Confirming Ion List | Select the confirming ions to be viewed. |
| Peak Labels | Displays or hides the peak labels (Label Area, Label Retention Time, Label Height, or Label Signal to Noise). |
| Show Peak Info | Displays peak information for the selected compound. For example: |
| | methyltestosterone |
| | Quan ion m/z: 267.10,285.20 Integration mode: Method |
| | Left RT: 1.87 Area: 684395 Apex RT: 1.99 Height: 95311 Right RT: 2.28 Noise: 530.01 Data file: steroids02 Filter: + c Full ms2 303.3@cid Detector: MS Trace: Mass range |
| | |
| Reset Scaling | Resets the original scaling after a zoom operation. |
| Peak Detection Settings | Opens the Peak Detection Settings dialog box for the selected compound. |

Qualification

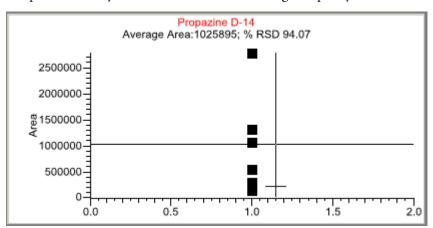
The Qualification pane displays the compound selected in the Quantification Data pane and Compounds pane.

The Qualification pane has five pages:

- Calibration Curve
- Spectra
- QED Spectra
- Confirming Ions
- Ion Overlay

Calibration Curve

The Calibration Curve page displays a graphical view of the calibration curve for the selected compound and key statistical values for evaluating the quality of the calibration.



Use the following procedures:

- To manually exclude a calibration point
- To zoom in on an area

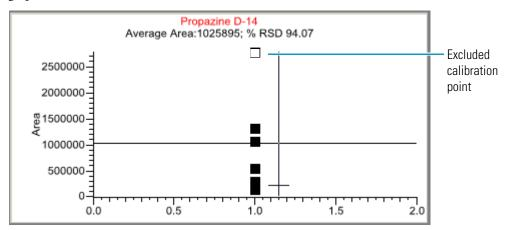
For a description of commands on the shortcut menu, see "Calibration Curve Shortcut Menu" on page 306.

❖ To manually exclude a calibration point

In the samples list, select the **Excluded** check box for the sample.

Use the horizontal scroll bar at the bottom of the table to scroll to the Excluded column.

When a value is no longer used for calibration, it is displayed as an empty box in the graphical view of the calibration curve.



❖ To zoom in on an area

- In the Calibration Curve plot, drag your cursor to delineate a rectangle around an area.
 The delineated area expands to fill the view.
- 2. To restore the method default view, right-click the calibration curve plot and choose **Reset Scaling** from the shortcut menu.

Changes to the calibration settings immediately apply to the entire results set, but you save them only when you save the batch.

Calibration Curve Shortcut Menu

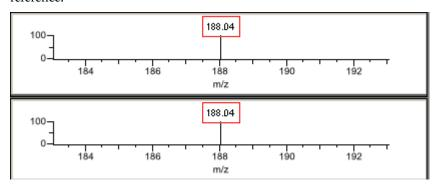
The Calibration Curve shortcut menu includes the following commands:

Table 65. Calibration Curve 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: Linear: Allows all settings with this exception: When Origin is set to Include, all Weighting values are grayed out and Weighting is set to Equal. Quadratic: Allows 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: Allows 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 | |
| Reset Scaling | Resets the original scale in the calibration curve pane. |

Spectra

The Spectra page displays a comparison of the spectra found in the data and the method reference.

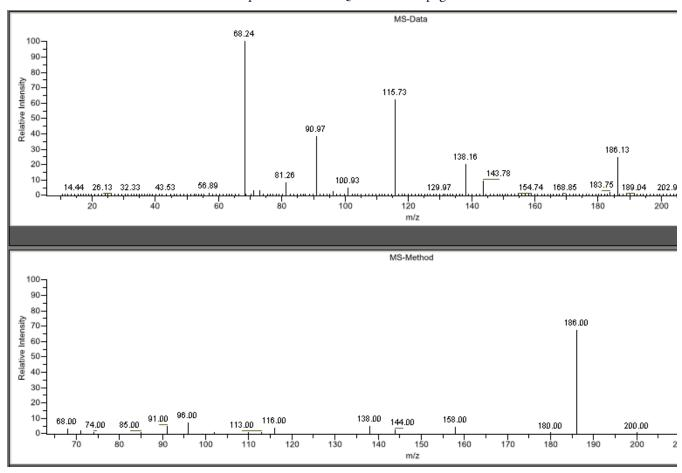


❖ To zoom in on an area

- 1. Drag your cursor to delineate a rectangle around an area.
 - The delineated area expands to fill the view.
- 2. To restore the method default view, right-click the spectra plot and choose **Reset Scaling** from the shortcut menu.

QED Spectra

The QED spectra display the averaged QED spectra from the raw data file and the datastore match. If the sample contains no QED data, the page is blank.



❖ To zoom in on an area

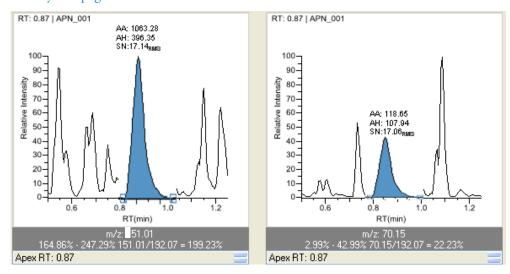
1. Drag your cursor to delineate a rectangle around an area.

The delineated area expands to fill the view.

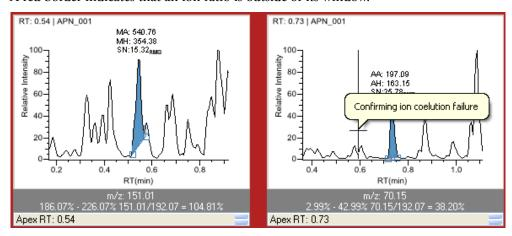
2. To restore the method default view, right-click the spectra plot and choose **Reset Scaling** from the shortcut menu.

Confirming Ions

The Confirming Ions page displays a graphical view of all qualifying/confirming ions for the selected sample and compound and displays calculated ion ratios and ion ratio acceptance windows. A red border indicates that an ion ratio is outside of its window. Depending on the method option settings, another ion view is available on the Ion Overlay page. See "Ion Overlay" on page 312.



A red border indicates that an ion ratio is outside of its window.



Use the following procedures:

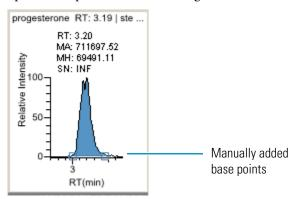
- To manually add a peak
- To remove a manually created peak
- To zoom in on a peak
- To change the displayed information for detected peaks
- To display the peak information window

For a description of commands on the shortcut menu, see "Confirming Ions Shortcut Menu" on page 311.

To manually add a peak

- Right-click the chromatogram plot and select **Add Peak** from the shortcut menu.
 If a peak is already detected, the Add Peak command is not enabled.
- 2. Click to indicate the left and right base points for the peak.

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 remove a manually created peak

Right-click the chromatogram plot, and choose **Remove Peak** from the shortcut menu. The application removes the manually added peak.

❖ To zoom in on a peak

- In the chromatogram plot, drag your cursor to delineate a rectangle around the peak.
 The delineated area expands to fill the view.
- 2. To restore the method default view, right-click the chromatogram plot and choose **Reset Scaling** from the shortcut menu.

To change the displayed information for detected peaks

- 1. Right-click the chromatogram plot and hold the cursor over **Peak Labels**.
- 2. Choose to display labels for the peak retention time, peak height, peak area, or signal to noise.
 - The label types in the list are selected for displayed labels and cleared for labels that are not displayed.
- 3. To remove a label, select the label type again and clear it.
 - Label settings are globally applied to quantification peaks, confirming peaks, and internal standard peaks.

Tip The labels do not always update on all peak displays. To update all labels, select a different compound, then reselect the compound whose labels you changed.

6 Using the Data Review Mode

Working in Data Review View

To display the peak information window

- 1. Right-click the quantification and confirming ion chromatogram plots.
- 2. Choose **Show Peak Info** from the shortcut menu.

Information in this window includes the data stream processed for this ion, consisting of a particular detector, filter (if applicable), and trace. For example, an MS detector might show a filter of "+ c Full ms [35.00-500.00]" and a Mass Range trace.

This window also shows the peak area and height, along with the retention times of the peak integration and the peak apex.

Confirming Ions Shortcut Menu

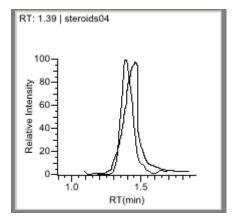
The Confirming Ions shortcut menu includes the following commands:

Table 66. Confirming lons shortcut menu commands

| Command | Description |
|---|---|
| Method Integration Settings | Displays method integration settings. |
| Manual Integration Settings | Displays manual integration settings. |
| Add Peak - Remove Peak - Cancel Add Peak | Adds a peak, removes a manually added peak, or cancels an add peak operation in progress. |
| Range Calc Method: Manual | Selects the method used to calculate the ion ratio range windows: Manual, Average, Weighted average, or Level. |
| Range Calc Level | Range based on the calibration level. |
| Target Ratio: | Specifies the theoretical ratio of the confirming ion's response to the quantification ion's response. |
| Window Type: | Specifies the Absolute or Relative calculation approach for determining the acceptable ion ratio range. |
| Window: % | Specifies the acceptable ion ratio range as a percentage. |
| Peak Labels | Displays or hides the peak labels (Label area, Label retention time, Label height, or Label signal to noise). |
| Show Peak Info | Displays peak information for the selected compound. For example: |
| | methyltestosterone |
| | Quan ion m/z: 267.10,285.20 Integration mode: Method |
| | Left RT: 1.87 Area: 684395 Apex RT: 1.99 Height: 95311 Right RT: 2.28 Noise: 530.01 Data file: steroids02 Filter: + c Full ms2 303.3@cid Detector: MS Trace: Mass range |
| Reset Scaling | Resets the original scaling after a zoom operation. |
| Peak Detection Settings | Opens the Peak Detection Settings dialog box for the selected compound. See "Peak Detection Settings dialog box" on page 302. |

Ion Overlay

The Ion Overlay page represents an overlay of the entire ion set—quantification and qualifying/confirming—for the selected sample and compound. Use this page to graphically review the peak apex alignment and co-eluting peak profiles.

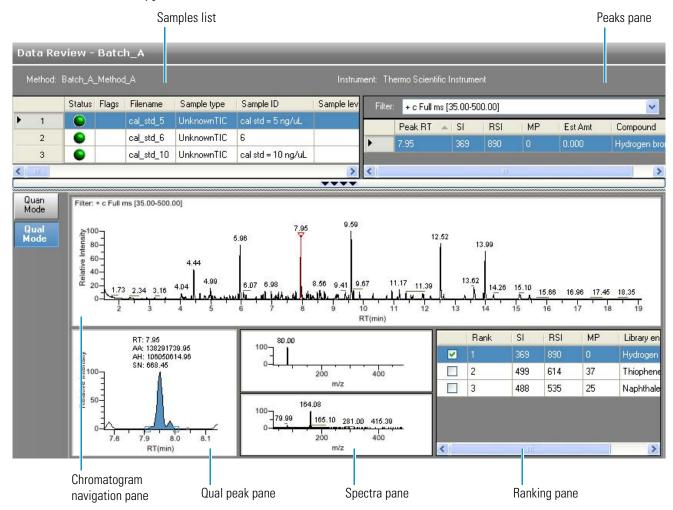


❖ To zoom in on a peak

- 1. Drag your cursor to delineate a rectangle around the peak.
 - The delineated area expands to fill the view.
- 2. To restore the method default view, right-click the chromatogram plot and choose **Reset Scaling** from the shortcut menu.

Qual Mode

Use the Qual Mode and the associated peaks pane to view qualitative information that complements the textual information for the selected Unknown/TIC sample. The Qual Mode view displays detected peaks for the selected sample. From the Qual Mode view, you can manually add peaks. The Qual Mode view is available only for Unknown/TIC sample types.



Tip To resize the panes, drag the separators that divide the panes.

In addition to the Samples list, the Qual Mode view displays data in the following panes:

- Peaks Pane
- Chromatogram Navigation Pane
- Qual Peak Pane
- Spectra Pane (Reference and Selected)
- Ranking Pane

Working in Data Review View

Peaks Pane

The peaks pane works with the samples list to display graphical values for a unique sample and peak combination. For detailed descriptions parameters on the peaks pane, see "Peaks pane" on page 316.

❖ To display peaks for a specific compound

1. From the samples list, select a sample.

Note If you choose a sample other than an Unknown/TIC sample, the TraceFinder application returns you to Quan Mode.

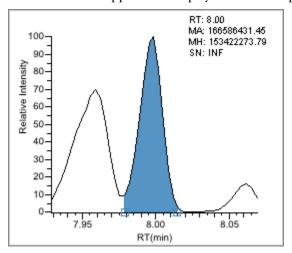
The peaks pane lists the retention times for peaks identified in the selected sample, the values for the best match methods for each peak, and the compound match.

The number of peaks that are listed is specified in the method. You can change the number of identified peaks in the Method Template Editor. See "Creating a Method Template" on page 188.

2. From the peaks pane, select a peak in the sample.

| | Peak RT (min) | SI | RSI | MP | Library entry |
|----------|---------------|-----|-----|----|--------------------------------------|
| | 5.01 | 844 | 847 | 33 | o-Toluidine |
| | 5.46 | 892 | 894 | 98 | 2-Cyclohexen-1-one, 3,5,5-trimethyl- |
|) | 8.00 | 937 | 937 | 80 | Acenaphthene |
| | 8.84 | 259 | 549 | 0 | Naphthalene, 1-methyl- |
| | 10.94 | 942 | 943 | 53 | Fluoranthene |

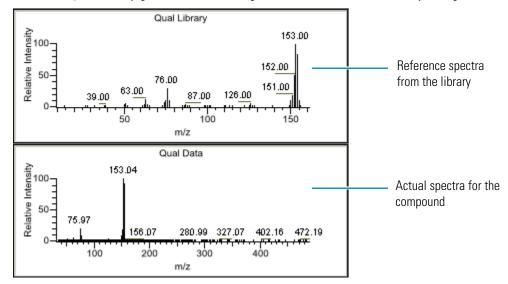
The TraceFinder application displays the selected peak in the qual peak pane.



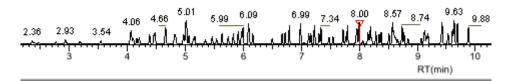
When you select a data-dependent sample, the peak can be from either a full scan or a QED spectrum of an SRM-filtered chromatogram.

The TraceFinder application displays the Spectra pane with two sections:

- The Qual Data pane shows spectra data for the peak in the raw data file.
- The Qual Library pane shows actual spectra for the identified library compound.



The TraceFinder application locates the selected peak in the navigation chromatogram.



❖ To remove a peak

- 1. Select a peak in the peaks pane.
- 2. Right-click and choose Remove Selected Peak from the shortcut menu.

The TraceFinder application removes the selected peak from the peaks list.

Note There is no undo for this action, but you can manually add a peak to redefine a removed peak. See "Chromatogram Navigation Pane" on page 317.

Working in Data Review View

Figure 77. Peaks pane

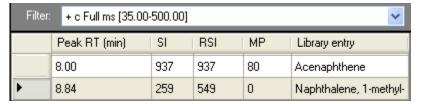


Table 67. Peaks pane parameters

| Command | 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": Filter: + c d Full ms2 179.15@cid35.00 [35.00-370.00] V Data-dependent filter | |
| Peak RT (min) | 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 method used to search the NIST library. | |
| MP | Match probability. | |
| Library entry | Library compound that matches the identified peak. | |
| Remove selected peak Shortcut menu command that removes the selected from the peaks list. | | |

Chromatogram Navigation Pane

The chromatogram navigation pane displays all peaks in the selected sample. The peak selected in the peaks pane displays a red marker.

For a description of commands on the shortcut menu, see "Chromatogram navigation pane shortcut menu commands" on page 318.

❖ To zoom in on a peak

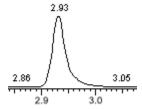
1. In the chromatogram navigation pane, drag the cursor to delineate a rectangle around the peak.

The delineated area expands to fill the view.

2. To restore the default view, right-click the chromatogram navigation pane and choose **Reset Scaling** from the shortcut menu.

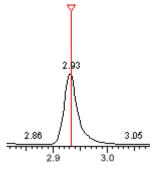
To manually add a peak

1. Zoom in to make it easier to identify the peak you want to add to the results set.



- 2. Right-click the chromatogram navigation pane, and choose **Add Peak** from the shortcut menu.
- 3. Click to indicate the left and right base points for the peak.

The TraceFinder application marks the peak in the chromatogram navigation pane.



The TraceFinder application places the peak delimiter tags at the base point locations and automatically updates the peak values in the peaks pane and qual peak pane.

RT: 2.93
MA: 53516847.82
MH: 29804603.78
SN: INF

Manually added base points

RT(min)

Figure 78. Qual peak pane with a manually added peak

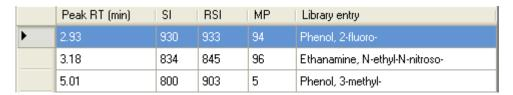


Figure 79. Chromatogram navigation pane

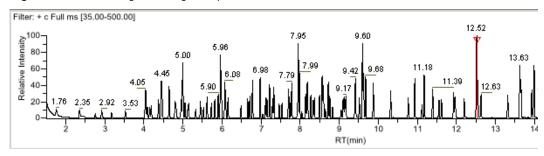


Table 68. Chromatogram navigation pane shortcut menu commands

| Command | Description |
|---------------|---|
| Add peak | Adds a peak. |
| Reset scaling | Resets the original scaling after a zoom operation. |

Qual Peak Pane

The qual peak pane displays the selected peak.

Use the following procedures:

- To zoom in on a peak
- To manually add a peak
- To remove a peak
- To switch between method and manual integration modes
- To change the displayed information for detected peaks

For a description of commands on the shortcut menu, see "Qual peak pane shortcut menu commands" on page 321.

To zoom in on a peak

- 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** from the shortcut menu.

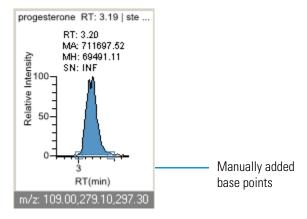
To manually add a peak

1. Right-click anywhere in the qual peak pane, and choose **Add Peak** from the shortcut menu.

If a peak is already detected, the Add Peak command is not enabled.

2. Click to indicate the left and right base points for the peak.

The TraceFinder 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 remove a peak

Right-click the chromatogram plot, and choose Remove Peak from the shortcut menu.

The TraceFinder application removes the peak displayed in the qual peak pane. All data for this peak are removed from the Qual Mode panes.

To switch between method and manual integration modes

Right-click the chromatogram view and choose **Method Integration** or **Manual Integration** from the shortcut menu.

Initially, the method 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, the 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 these data identify the manual modifications.

❖ To change the displayed information for detected peaks

- 1. Right-click the chromatogram plot and hold the cursor over **Peak Labels**.
- 2. Choose to display labels for the peak retention time, peak height, peak area, or signal to noise.

The label types in the list are selected for displayed labels and cleared for labels that are not displayed.

3. To remove a label, select the label type again and clear it.

Label settings are globally applied to qual peaks, confirming peaks, and internal standard peaks.

Tip The labels do not always update on all peak displays. To update all labels, select a different compound, and then reselect the compound whose labels you changed.

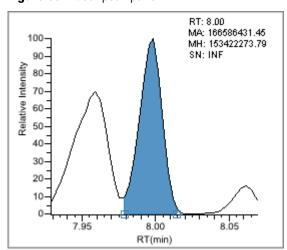


Figure 80. Qual peak pane

The qual peak pane shortcut menu includes the following commands:

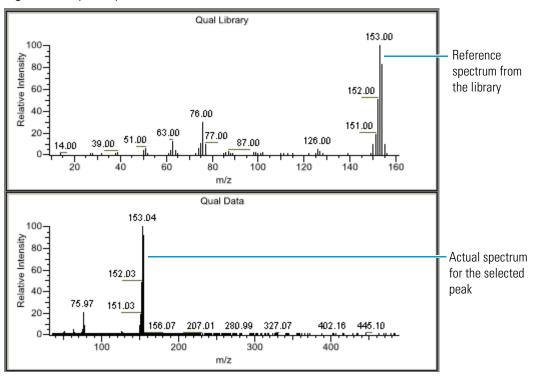
Table 69. Qual peak pane shortcut menu commands

| 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 Signal to Noise). |
| Remove peak | Removes the peak displayed in the Qual pane. |

Spectra Pane (Reference and Selected)

The Spectra pane displays the reference spectra and the spectra for the selected sample. The top pane displays the reference spectra for the identified compound from the library; the bottom pane displays the spectra for the selected peak.

Figure 81. Spectra pane



To zoom in on a peak

- In the spectra 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** from the shortcut menu.

Acenaphthene

Ranking Pane

The ranking pane displays the three 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:

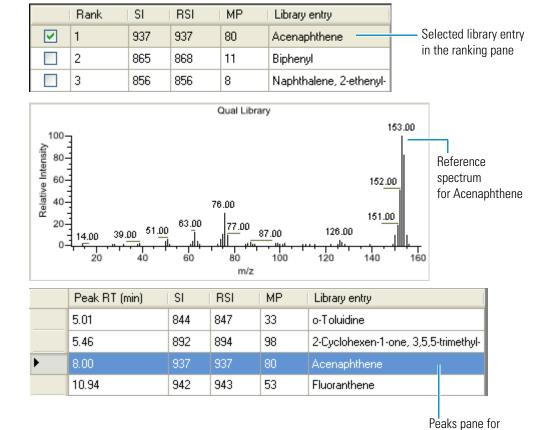


For a detailed description of ranking pane parameters, see "Ranking pane" on page 324.

❖ To change the library entry for a selected peak

In the ranking pane, select the check box for the library entry you want to use to identify the selected peak.

- In the Spectra pane, the reference spectra change to show the spectra for the selected library entry.
- In the peaks pane, the SI, RSI, MP, and Library Entry values update to reflect the selected library entry.



6 Using the Data Review Mode

Working in Data Review View

Figure 82. Ranking pane

| | Rank | SI | RSI | MP | Library entry |
|---|------|-----|-----|----|-------------------------|
| ~ | 1 | 937 | 937 | 80 | Acenaphthene |
| | 2 | 865 | 868 | 11 | Biphenyl |
| | 3 | 856 | 856 | 8 | Naphthalene, 2-ethenyl- |

Table 70. Ranking pane parameters

| Command | Description |
|----------------------------------|--|
| <check box="" column=""></check> | 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 method used to search the NIST library. |
| RSI | Reverse search index method used to search the NIST library. |
| MP | Match probability. |
| Library entry | Library compound that matches the identified peak. |

Working in Report View

Use the Report View to display or generate reports for the currently selected batch in the Data Review mode. You must process each sample in the batch before you can view or generate a sample-level report for that sample.

This section includes the following topics:

- Viewing Reports
- Generating Reports
- Working with Reports
- Working with Active View

Figure 83. Report View in Data Review mode



• View Only: Displays a PDF or Excel spreadsheet preview of the selected report type for the batch, sample, or compound. See "Viewing Reports" on page 326.

Preview reports for all Standard report types are always available. You must generate Custom and Target Screening report types before they are available in this list.

The Report View page displays one of the following report outputs:

- Standard reports as PDF files
- Custom reports in XLSM format
- Target Screening reports as PDF files
- Generate Only: Creates all specified report output formats for the selected sample- or batch-level report. See "Generating Reports" on page 330.

❖ To open the Report View

- 1. Click **Data Review** in the navigation pane from any mode.
- 2. In the Data Review navigation pane, click **Report View**.



The Report View for the currently selected batch opens.

Viewing Reports

Use the View Only features to view all configured standard reports and any custom or target screening reports that you have generated. After you generate a report, the application displays the report in the View Only report list.

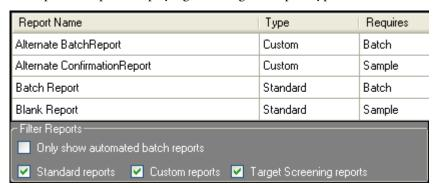
Use the following procedures:

- To select a report
- To select a sample
- To select a compound
- To select a sample and a compound

❖ To select a report

- 1. Select the **View Only** option.
- 2. Click the Select a Report box.

The report list opens, displaying all configured report types.



These reports reflect the Displayed Reports selections in the Configuration mode. To change the configured reports that are available in this view, see "Specifying the Reports Configuration" on page 58.

To sort the reports, click the column headers. The application maintains this sort order each time you open the Report View for this batch.

To help organize your reports, you can filter the list.

3. To limit the types of reports you want to display in the report list, select any combination of report filter options in the Filter Reports area.

Table 71. Filter Reports options

| Option | Behavior |
|-----------------------------------|--|
| Only show automated batch reports | Displays only reports that have an output format specified in the Automated Batch Reports area in the Batch View. See "Editing Report Output Formats" on page 274. |
| Standard reports | Displays Standard report types. |
| Custom reports | Displays all generated Custom report types. Custom reports are not available for viewing until you have generated the report. |
| Target Screening reports | Displays all generated Target Screening reports. Target Screening reports are not available for viewing until you have generated the report. |

Note When you make changes to the method in the Local Method view, to the peaks in the Data Review view, or to the samples in the Batch View, you must regenerate the custom or target screening reports to see the effects of those changes.

4. Double-click the name of the report.

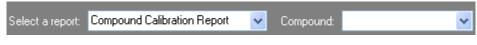
The report list closes.

- When the selected report is a batch-level report, the application displays the report on the Report View page.
- When the selected report includes separate reports for each sample, you must select a sample file.



Follow the procedure "To select a sample" on page 328.

• When the selected report includes separate reports for each compound, you must select a compound.



Follow the procedure "To select a compound" on page 328.

• When the selected report includes separate reports for each sample and each compound in the sample, you must select both a sample and a compound.

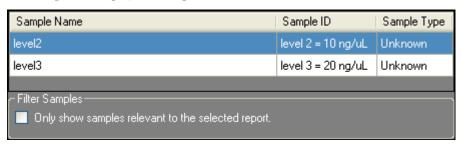


Follow the procedure "To select a sample and a compound" on page 329.

❖ To select a sample

1. Click the Sample File box.

The sample list displays all samples in the batch.



2. To show only samples that would be included in the selected report, select the **Only Show Samples Relevant...** check box.

For example, if you selected the Check Standard Report, the sample list displays only Chk Std samples.

Note Click the column headers to sort the samples. The application maintains this sort order each time you open the Report View for this batch.

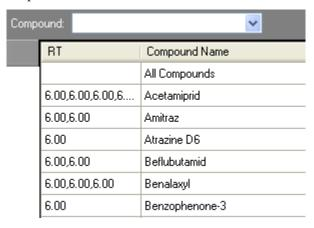
3. Double-click the name of the sample.

The sample list closes. The Report View page displays the sample-level report.

To select a compound

1. Click the Compound list.

The compound list displays the names and retention times of all compounds in the sample.



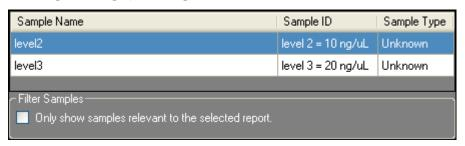
2. Double-click a single compound or **All Compounds**.

The compound list closes. The Report View page displays the compound-level report.

❖ To select a sample and a compound

1. Click the Sample File box.

The sample list displays all samples in the batch.



2. To show only samples that would be included in the selected report, select the **Only Show Samples Relevant...** check box.

For example, if you selected the Check Standard Report, the sample list displays only Chk Std samples.

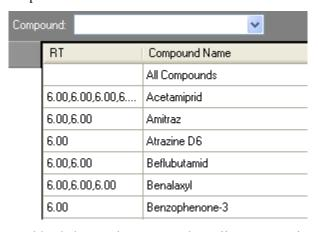
Tip Click the column headers to sort the samples. The application saves this sort order in the Report View for this batch.

3. Double-click the name of the sample.

The sample list closes.

4. Click the Compound list.

The compound list displays the names and retention times of all compounds in the sample.



5. Double-click a single compound or All Compounds.

The compound list closes.

The Report View page displays the compound-level report for the selected sample and compound.

Generating Reports

Use the Generate Only features to create sample-level reports. You cannot use the View Only features to view custom or target screening reports until you generate the report. When you make changes to the method in the Local Method view or to the peaks in the Data Review view, you must regenerate the custom or target screening reports to see the effects of those changes.

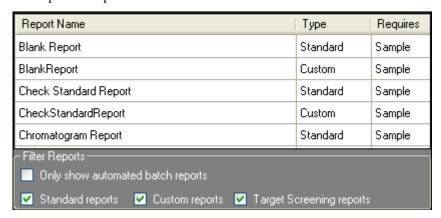
Use the following procedures:

- To select a report
- To select a sample

❖ To select a report

- 1. Select the **Generate Only** option.
- 2. Click the Select a Report box.

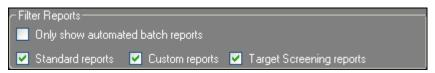
The report list opens.



All configured sample-level report types are displayed in the list. You cannot generate batch-level or compound-level reports from this view. To change the configured reports that are available in this view, see "Specifying the Reports Configuration" on page 58.

If you have many reports, you can filter the list.

3. To limit the types of reports you want to display in the report list, select any combination of report filter check boxes in the Filter Reports area.

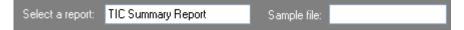


| Option | Behavior |
|-----------------------------------|---|
| Only show automated batch reports | Displays only sample-level reports that have an output format specified in the Automated Batch Reports area in the Batch View. See "Editing Report Output Formats" on page 274. |
| | If you have only batch-level reports specified in the Batch View, selecting this option excludes all reports in the Report Name list. |
| Standard reports | Displays sample-level Standard report types. |
| Custom reports | Displays sample-level Custom report types. |
| Target Screening reports | Displays sample-level Target Screening report types. |

Note Click the column headers to sort the samples. The application saves this sort order in the Report View for this batch.

4. Double-click the name of the report.

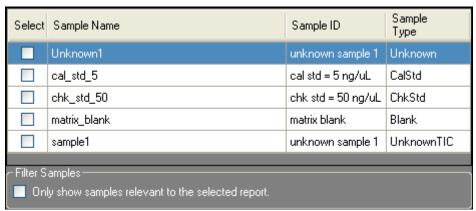
The report list closes. You must select a sample file for the selected report.



❖ To select a sample

1. Click the Sample File box.

The sample list displays all samples in the batch.



2. To show only samples that would be included in the selected report, select the **Only Show Samples Relevant...** check box.

6 Using the Data Review Mode

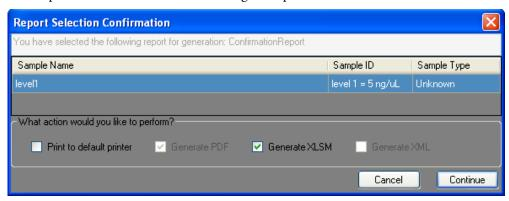
Working in Report View

For example, if you selected the Check Standard Report, the sample list displays only Chk Std samples.

Note Click the column headers to sort the samples. The application saves this sort order in the Report View for this batch.

- 3. Select the check box for each sample you want to include in the report.
- 4. Click Generate.

The Report Selection Confirmation dialog box opens.

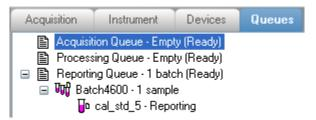


5. In the What Action Would You Like to Perform area, select the types of reports you want to create.

Note The application automatically selects required output formats. These options are not editable.

6. Click Continue.

The application submits the selected samples to the report queue.



When you have already generated this report in the Batch View or Acquisition Mode, the new report is time-stamped to differentiate it from the original report.

7. To view the report you generated, follow the instructions in "Viewing Reports" on page 326.

Note When you make changes to the method in the Local Method view, to the peaks in the Data Review view, or to the samples in the Batch View, you must regenerate the custom or target screening reports to see the effects of those changes.

Working with Reports

Use the buttons on the Report View page to view, print, or export a report.

- A PDF report view is available for all Standard and Target Screening report types.
- An Excel Macro-Enabled Workbook report view is available for any Custom report types
 that you have generated with the Generate XLSM option selected.

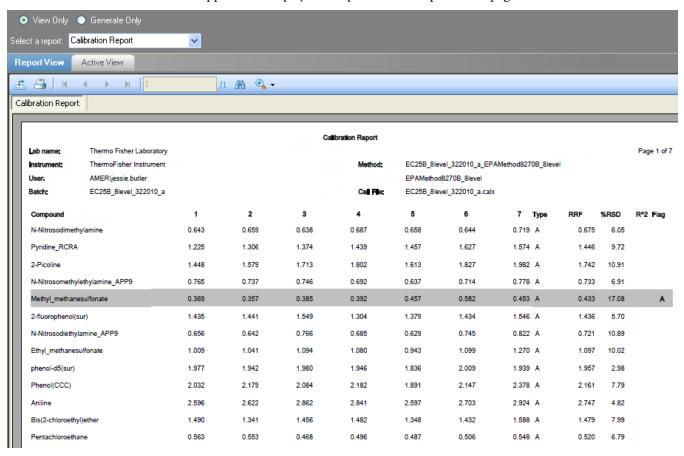
Use the following procedures:

- To print a standard or target screening report
- To export a standard report
- To search for text
- To enlarge the report text

To print a standard or target screening report

- 1. Select the report you want to print from the Select a Report list.
- 2. (Optional) Select a sample from the Sample File list.

The application displays the report on the Report View page.



6 Using the Data Review Mode

Working in Report View

3. Click the **Print Report** button,

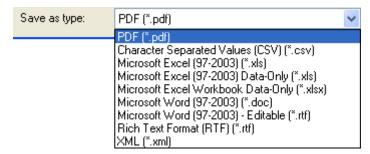


The Print dialog box for your default printer opens.

4. Follow the typical procedure to print from your printer. Landscape reports automatically rotate to fit the paper.

To export a standard report

- 1. Select the report you want to print from the Select a Report list.
- 2. (Optional) Select a sample from the Sample File list. The application displays the report on the Report View page.
- 3. Click the **Export Report** button, <u></u> The Export Report dialog box opens.
- 4. Locate the folder where you want to write the report file.
- 5. Type a file name for the exported report file.
- 6. Select a file type from the Save as Type list:



7. Click Save.

The TraceFinder application saves the file as the specified file type and writes the report file to the specified folder.

To search for text

- 1. Select a report from the Select a Report list.
- 2. (Optional) Select a sample from the Sample File list.

The application displays the report on the Report View page.



The Find Text dialog box opens.

4. Enter your text and click **Find Next**.

When the TraceFinder application locates the text, it encloses the text in a red box.

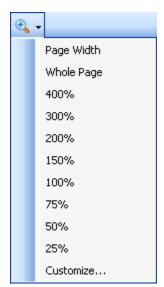
Sample ID APN001 APN002 APN003

To enlarge the report text

- 1. Select a report from the Select a Report list.
- 2. (Optional) Select a sample from the Sample File list.

The application displays the report on the Report View page.

3. Click the **Zoom** button, , and select a zoom scale.



Working with Active View

Use the Active View page to view quantitative data for each sample in a report. Data in the Active View are labeled with flag information. These flags are based on a comparison of the batch data to criteria defined in the master method.

❖ To display the Active View page

Click the **Active View** tab.

The Active View page displays quantitative data and QAQC error flags for each sample. See "Active View page" on page 337.

To display a report

1. Select a report type from the Select a Report list.

Only the report types created for the current batch are displayed in the list.

2. (Optional) When the report type includes separate reports for each sample, select a sample file.



❖ To filter which compounds to display

Click the Showing button to display either all compounds or only compounds that are flagged for failing a QAQC test.



Figure 84. Active View page

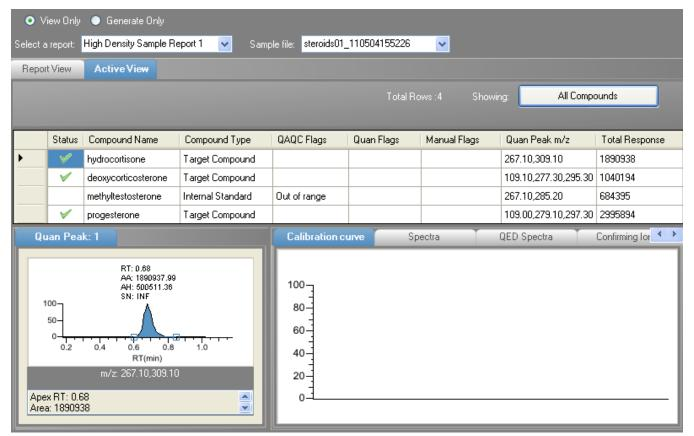


Table 72. Active View parameters (Sheet 1 of 4)

| Parameter | Description |
|-----------------|--|
| View Only | The Active View pane is available only when View Only is selected. |
| Select a report | Displays the report types created for the current batch. |
| Sample file | Used when the report type includes separate reports for each sample. |
| Total Rows | The number of compound rows currently displayed in the pane. |
| Showing | Displays all compounds or only the flagged compounds. |

6 Using the Data Review Mode Working in Report View

Table 72. Active View parameters (Sheet 2 of 4)

| Parameter | Description |
|-----------------------|--|
| Column headings | Many column headings are specific to individual reports. See "Active View Report Contents" on page 341. |
| Status | Indicates the status of the reported compound. |
| | A yellow check mark indicates one of the following conditions: |
| | The compound was manually integrated. |
| | Any of the confirming peaks was manually integrated. |
| | - The compound has quan flags. |
| | A red check mark indicates that the QAQC checks failed. |
| | A green check mark indicates that none of these conditions exist. |
| | When the compound is an internal standard, warnings are displayed only on the internal standard report. The Status column is blank for Manual Integration reports. |
| Compound Name | Alphanumeric name assigned to the compound. |
| Compound Type | Target Compound, Internal Standard, or Surrogate. |
| QAQC Flags | Indicates that the QAQC check for the sample failed. |
| | The QAQC column is not used for Manual Integration reports. |
| Quan Flags | • Limit of Detection (LOD) |
| | Limit of Quantitation (LOQ)Limit of Reporting (LOR) |
| | Values between the limit of detection and the limit of quantitation, known as the J flag Upper Limit of Linearity (ULOL) |
| | Quan flags do not apply to these sample types: Cal Std, Chk Std, Matrix Blank, or Solvent. |
| | The Quan Flag column is not used for Manual Integration reports. |
| Manual Flags | Indicates manually integrated peaks. |
| | • M indicates a manually integrated quan peak. |
| Depending on the sele | m indicates a manually integrated confirming peak. cted report, the Active View page contains any or all of the following parameters: |
| Quan Peak m/z | Mass-to-charge ratio for the selected quantitation peak. |
| Total Response | The sum of all Quan Peak Response values for the compound. |
| Quan Peak Response | Response of the quan peak. |
| Quan peak RT | Retention time for the quan peak. |
| Theoretical amount | Theoretical amount of the compound. Reports N/A when not applicable. |
| Concentration | |
| | Mass of the confirming peak. |

Table 72. Active View parameters (Sheet 3 of 4)

| Parameter | Description |
|---------------------------------------|--|
| Confirming <i>n</i> Response | Response of the confirming peak. |
| Confirming <i>n</i> Manual Flag | Indicates a manually integrated confirming peak. |
| Confirming <i>n</i> Ion Ratio Flag | Indicates that the ion ratio is out of range. |
| Confirming <i>n</i> Ion Ratio | Actual ratio of the confirming ion response to the quan ion response. |
| Confirming n Range | Acceptable range for the confirming ion. |
| Retention Time | The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| Quan mass | The mass-to-charge ratio used to determine the peak area and peak height of the compound. |
| Response | Sum of all Quan Peak Response values for the compound. |
| Injection concentration | Calculated amount as the sample was injected, with no conversion applied. |
| Injection Units | Injection units specified on the Calibration page in Method Development mode. See "Calibration" on page 151. |
| Sample Concentration | The injected concentration multiplied by the conversion factor. |
| Sample Units | Sample units specified on the Calibration page in Method Development mode. See "Calibration" on page 151. |
| QIon | Mass range for the quan peak. |
| RT | Retention time. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| Manual Integration repo | rts |
| m/z | Mass-to-charge ratio for the quan peak. |
| Method RT | Apex retention time for the method-integrated peak. |
| Method Peak Height | Height of the method-integrated peak. |
| Method Peak Area | Area of the method-integrated peak. |
| Manual RT | Apex retention time for the manually integrated peak. |
| Manual Peak Height | Height of the manually integrated peak. |
| Manual Peak Area | Area of the manually integrated peak. |

6 Using the Data Review Mode Working in Report View

Table 72. Active View parameters (Sheet 4 of 4)

| Parameter | Description | |
|---------------------------|--|--|
| Internal Standard reports | | |
| Std Response | Average of the internal standard's response as found in the calibration file. | |
| Minimum Response | Minimum response time as specified on the ISTD page in Method Development mode. See "ISTD" on page 166. | |
| Maximum Response | Maximum response time as specified on the ISTD page in Method Development mode. See "ISTD" on page 166. | |
| Sample Response | Area found in the sample. | |
| Std RT | Average retention time as found in the calibration file. | |
| Min RT | Minimum retention time as specified on the ISTD page in Method Development mode. See "ISTD" on page 166. | |
| Max RT | Maximum retention time as specified on the ISTD page in Method Development mode. See "ISTD" on page 166. | |
| Sample RT | Retention time found in the sample. | |
| Graphical data | | |
| Quan Peak 1 | | |
| Calibration curve | Displays a graphical view of the calibration curve for the selected compound and key statistical values for evaluating the quality of the calibration. | |
| Spectra | Displays a comparison of the spectra found in the data and the method reference. | |
| QED Spectra | Displays the averaged QED spectra from the raw data file and the datastore match. If the sample contains no QED data, the page is blank. | |
| Confirming Ions | Displays a graphical view of all qualifying/confirming ions for the selected sample and compound, and displays calculated ion ratios and ion ratio acceptance windows. | |

Active View Report Contents

Each standard report that uses the Active View displays values that are common to all reports:

• Common Active View report columns

In addition to the common values, the following reports display additional active view features:

- Blank Report Active View columns
- Calibration Report Active View columns
- High Density Sample Report 1 and High Density Sample Report 1 Long Active View columns
- High Density Sample Report 2 and High Density Sample Report 2 Long Active View columns
- High Density Sample Report 3 and High Density Sample Report 3 Long Active View columns
- Internal Standard Summary Report Active View columns
- Ion Ratio Failure Report Active View columns
- Manual Integration Report Active View columns
- LCSLCSD Report Active View columns (Environmental and Food Safety only)
- Method Detection Limit Report Active View columns (Environmental and Food Safety only)
- Method Validation Report Active View columns (Environmental and Food Safety only)
- MSMSD Report Active View columns (Environmental and Food Safety only)
- Quantitation Report Active View columns
- Solvent Blank Report Active View columns

Table 73. Common Active View report columns

| Column | Description |
|---------------|--|
| Status | Indicates the status of the reported compound. |
| | A yellow caution sign indicates one of the following conditions: |
| | The compound was manually integrated. |
| | Any of the confirming peaks was manually integrated. |
| | The compound has quan flags. |
| | The compound has a QAQC failure. |
| | A green check mark indicates that none of these conditions exists. |
| | When the compound is an internal standard, warning flags are displayed only on the internal standard report. |
| Compound name | Alphanumeric name assigned to the compound. |
| Compound type | Target Compound, Internal Standard, or Surrogate. |
| QAQC flags | Indicates that the QAQC check for the sample failed. |
| | This column is not included on the Method Validation and MDL reports. |
| Quan flags | Limit of Detection (LOD) Limit of Quantitation (LOQ) Limit of Reporting (LOR) Values between the limit of detection and the limit of quantitation, known as the J flag Upper Limit of Linearity (ULOL) |
| | Quan flags do not apply to these sample types: Cal Std, Chk Std. Matrix Blank, or Solvent. |
| | This column is not included on the Calibration report. |
| | This column is not included on the Calibration Curve report. |
| | This column is not included on the Method Validation, Method Detection Limit, LCS, or LCSD reports. |
| Manual flags | Indicates manually integrated peaks. M indicates a manually integrated quan peak. m indicates a manually integrated confirming peak. |

Table 74. Blank Report Active View columns

| Column | Description |
|----------------|--|
| Retention Time | Retention time for the quan mass. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| Quan Mass | Mass range for the quan peak. |
| Response | Sum of all Quan Peak Response values for the compound. |
| Inj Conc | Calculated amount as the sample was injected, with no conversion applied. |
| Inj Units | Injection units specified on the Calibration page in Method Development mode. See "Calibration" on page 151. |
| Sample Conc | Calculated amount multiplied by the conversion factor. |
| Sample Units | Sample units specified on the Calibration page in Method Development mode. See "Calibration" on page 151. |

Table 75. Calibration Report Active View columns (Sheet 1 of 2)

| Column | Description |
|------------------|---|
| Curve Type | The type of curve used when calibrating the compound (linear, quadratic, or average response factor). |
| Average RF | The average response factor. Applicable if curve type is Average RF. |
| Average Response | The average response for the internal standard across all calibration points. Applies only to Internal Standard sample types. |
| A0 | The value with no X. Applies only to linear and quadratic curves. |
| A1 | The X value. Applies only to linear and quadratic curves. |
| A2 | The X^2 value. Applies only to quadratic curves. |
| R^2 | The minimum correlation coefficient (r^2) for an acceptable calibration (when in linear or quadratic mode). |

Table 75. Calibration Report Active View columns (Sheet 2 of 2)

| Column | Description |
|--------|---|
| RSD | Relative standard deviation. Applies only to internal standards and targets calibrated with an average RF curve. |
| Level | The column specifies the level name; the field value specifies the data point used in calibration. This field can be Response Factor for external calibration, Response Ratio for internal linear or quadratic, or Relative Response Factor for Internal Average RF. There is one column for each level in the curve. If the batch uses an extended calibration, there might be more columns than calibration standards in the current batch. |

Table 76. High Density Sample Report 1 and High Density Sample Report 1 Long Active View columns

| Column | Description |
|--------------------|--|
| m/z | Mass-to-charge ratio for the quan peak. |
| Total Response | The sum of all Quan Peak Response values for the compound. |
| Quan Peak Response | Response of the quan peak. |
| Quan Peak RT | Retention time for the quan peak. |
| T Amount | Theoretical amount of the compound. Reports N/A when not applicable. |
| Conc | Calculated (injected) amount. |

Table 77. High Density Sample Report 2 and High Density Sample Report 2 Long Active View columns (Sheet 1 of 2)

| Column | Description |
|--------------------------|--|
| m/z | Mass-to-charge ratio for the quan peak. |
| Total Response | Sum of all Quan Peak Response values for the compound. |
| Quan Peak Response | Response of the quan peak. |
| Quan Peak RT | Retention time for the quan peak. |
| T Amount | Theoretical amount of the compound. Reports N/A when not applicable. |
| Conc | Calculated (injected) amount. |
| Confirming 1 Mass | Mass of the confirming peak. |
| Confirming 1 Response | Response of the confirming peak. |
| Confirming 1 Manual Flag | Indicates a manually integrated confirming peak. |

Table 77. High Density Sample Report 2 and High Density Sample Report 2 Long Active View columns (Sheet 2 of 2)

| Column | Description |
|-----------------------------|---|
| Confirming 1 Ion Ratio Flag | Indicates that the ion ratio is out of range. |
| Confirming 1 Ion Ratio | Actual ratio of the confirming ion response to the quan ion response. |
| Confirming 1 Range | Acceptable range for the confirming ion. |

Table 78. High Density Sample Report 3 and High Density Sample Report 3 Long Active View columns

| Column | Description |
|-----------------------------|---|
| m/z | Mass-to-charge ratio for the quan peak. |
| Total Response | Sum of all Quan Peak Response values for the compound. |
| Quan Peak Response | Response of the quan peak. |
| Quan Peak RT | Retention time for the quan peak. |
| T Amount | Theoretical amount of the compound. Reports N/A when not applicable. |
| Conc | Calculated (injected) amount. |
| Confirming 1 Mass | Mass of the confirming peak. |
| Confirming 1 Response | Response of the confirming peak. |
| Confirming 1 Manual Flag | Indicates a manually integrated confirming peak. |
| Confirming 1 Ion Ratio Flag | Indicates that the ion ratio is out of range. |
| Confirming 1 Ion Ratio | Actual ratio of the confirming ion response to the quan ion response. |
| Confirming 1 Range | Acceptable range for the confirming ion. |
| Confirming 2 Mass | Mass of the confirming peak. |
| Confirming 2 Response | Response of the confirming peak. |
| Confirming 2 Manual Flag | Indicates a manually integrated confirming peak. |
| Confirming 2 Ion Ratio Flag | Indicates that the ion ratio is out of range. |
| Confirming 2 Ion Ratio | Actual ratio of the confirming ion response to the quan ion response. |
| Confirming 2 Range | Acceptable range for the confirming ion. |

Table 79. Internal Standard Summary Report Active View columns

| Column | Description |
|------------------|--|
| Std Response | Average of the internal standard's response as found in the calibration file. |
| Minimum Response | Minimum response time as specified on the ISTD page in Method Development mode. See "ISTD" on page 166. |
| Maximum Response | Maximum response time as specified on the ISTD page in Method Development mode. See "ISTD" on page 166. |
| Sample Response | Area found in the sample. |
| Std RT | Average retention time as found in the calibration file. |
| Min RT | Minimum retention time as specified on the ISTD page in Method Development mode. See "ISTD" on page 166. |
| Max RT | Maximum retention time as specified on the ISTD page in Method Development mode. See "ISTD" on page 166. |
| Sample RT | Retention time found in the sample. |

Table 80. Ion Ratio Failure Report Active View columns

| Column | Description |
|-------------------|--|
| Quan Ion | The ion for quan peak. |
| Qual Ion | The ion for the confirming peak. |
| Quan Ion Response | Response of the quan ion. |
| Qual Ion Response | Response of the qual ion. |
| Ratio | The ratio of the confirming ion response to the quan ion response. |
| Range | The acceptable range. |

Table 81. Manual Integration Report Active View columns

| Column | Description |
|--------------------|---|
| m/z | Mass-to-charge ratio for the quan peak. |
| Method RT | Apex retention time for the method-integrated peak. |
| Method Peak Height | Height for the method-integrated peak. |
| Method Peak Area | Area for the method-integrated peak. |
| Manual RT | Apex retention time for the manually integrated peak. |
| Manual Peak Height | Height of the manually integrated peak. |
| Manual Peak Area | Area of the manually integrated peak. |

Table 82. LCSLCSD Report Active View columns (Environmental and Food Safety only)

| Column | Description |
|------------------------|--|
| Spike Amount | Lab control theoretical concentration. |
| LCS Concentration | Lab control spike concentration. |
| LCS % Received | Lab control concentration percentage. |
| Lower Limit % | Recovery lower limit as specified in the master method. See the Limits section in "Editing the Flags Page" on page 159. |
| Upper Limit % | Recovery upper limit as specified in the master method. See the Limits section in "Editing the Flags Page" on page 159. |
| LCSD Concentration | Lab control spike duplicate concentration. |
| LCSD % Received | Lab control spike duplicate concentration percentage. |
| RPD | Lab control relative percentage difference. |
| Max RPD | Lab control spike maximum relative percentage difference (set in method). |
| Number of Rec Failures | Number of recovery failures for lab control spike concentration and lab control spike concentration duplicate. |
| Number of RPD Failures | Number of relative percentage difference failures for lab control spike concentration and lab control spike concentration duplicate. |

Note For LCSLCSD batch reports, the application displays the active view peak graphics only when you click a field pertaining to a sample, such as the LCS or LCSD concentration fields.

Table 83. Manual Integration Report Active View columns

| Column | Description |
|--------------------|---|
| m/z | Mass-to-charge ratio for the quan peak. |
| Method RT | Apex retention time for the method-integrated peak. |
| Method Peak Height | Height for the method-integrated peak. |
| Method Peak Area | Area for the method-integrated peak. |
| Manual RT | Apex retention time for the manually integrated peak. |
| Manual Peak Height | Height of the manually integrated peak. |
| Manual Peak Area | Area of the manually integrated peak. |

Table 84. Method Detection Limit Report Active View columns (Environmental and Food Safety only)

| Column | Description |
|----------|--|
| Avg Conc | The average of the concentration for the compound across all samples. |
| Std Dev | The standard deviation of the concentration. |
| t-stat | The t-statistic value defined as the ratio of a coefficient to its standard error. |
| % RSD | %RSD of concentrations |
| MDL | Method detection limits. The calculated limit of detection. |

Note For Method Detection Limit batch reports, the application displays the active view peak graphics only when you click a field pertaining to a sample. These numbered fields are to the right of the MDL column.

Table 85. Method Validation Report Active View columns (Environmental and Food Safety only)

| Column | Description |
|--|--|
| Avg Conc | The average of the concentration for the compound across all samples. |
| Theo Conc | Values for each compound that represent the expected theoretical concentration of that compound in the sample as defined in the master method. See the Meth Val section in "Editing the Flags Page" on page 159. |
| % Diff | The percentage difference calculated as ([MethodValidationMeanValue minus the Theo Conc] divided by the Theo Conc) multiplied by 100. |
| Min Conc | Calculated by (Min recovery percent * Theo Conc) divided by 100. |
| Max Conc | Calculated by (Max recovery percent * Theo Conc) divided by 100. |
| % RSD | %RSD of concentrations |
| Max % RSD | The maximum relative standard deviation (RSD) of the set of observed concentrations for a component across the set of method validation samples (when in average RF mode) as defined in the master method. See the Meth Val section in "Editing the Flags Page" on page 159. |
| Calculated Amount <i><sample< i=""> Name></sample<></i> | This field is reproduced for every Method Val sample. |

Table 86. MSMSD Report Active View columns (Environmental and Food Safety only)

| Column | Description |
|------------------------|--|
| Unknown Concentration | Concentration of the unknown sample. |
| Spike Amount | Matrix spike theoretical concentration. |
| MS Concentration | Matrix spike concentration. |
| MS % Received | Matrix spike concentration percentage. |
| Lower Limit % | Recovery lower limit as specified in the master method. See the Limits section in "Editing the Flags Page" on page 159. |
| Upper Limit % | Recovery upper limit as specified in the master method. See the Limits section in "Editing the Flags Page" on page 159. |
| MSD Concentration | Matrix spike duplicate concentration. |
| MSD % Received | Matrix spike duplicate concentration percentage. |
| RPD | Matrix spike relative percentage difference. |
| Max RPD | Maximum relative percentage difference as specified in the master method. See the Lab Control section in "Editing the Flags Page" on page 159. |
| Number of Rec Failures | Number of matrix spike and matrix spike duplicate failures. |
| Number of RPD Failures | Number of relative percentage difference failures. |

Note For MSMSD batch reports, the active view peak graphics are shown only when you click a field pertaining to a sample, such as Unknown, MS, or MSD concentration fields.

Table 87. Quantitation Report Active View columns (Sheet 1 of 2)

| Column | Description |
|----------|---|
| RT | Retention time for the peak. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| QIon | Mass range for the quan peak. |
| Response | Sum of all quan peak response values for the compound. |

Table 87. Quantitation Report Active View columns (Sheet 2 of 2)

| Column | Description |
|------------------------|---|
| Injected Concentration | Calculated amount as the sample was injected, with no conversion applied. |
| | As each additional sample is processed, calibration data change; therefore, except for the final sample in a batch, a report in active view or report view shows different values from a physical (PDF, XML, or printed) report created at the end of processing. To avoid this discrepancy, do one of the following: • For the standard Quantitation Report or Quantitation Report - 2, observe the active or report view for only the last sample in the batch. • For the custom Quantitation Report, make the report a batch-level report. |
| Injected units | Injection units specified on the Calibration page in Method Development mode. See "Calibration" on page 151. |
| Sample Conc | Calculated injection amount multiplied by the conversion factor. See the Injected Concentration description. |
| Sample units | Sample units specified on the Calibration page in Method Development mode. See "Calibration" on page 151. |

Table 88. Solvent Blank Report Active View columns

| Column | Description |
|-------------|--|
| RT | Retention time for the quan peak. The time after injection when the compound elutes. The total time that the compound is retained on the column. |
| QIon | Mass range for the quan peak. |
| Response | Sum of all Quan Peak Response values for the compound. |
| Method | Method of evaluation defined in the method. |
| Upper Limit | Defined in the method. |

Working in the Local Method View

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 *BatchName MasterMethodName*.

❖ To open the Local Method View

1. Do one of the following:

From the dashboard, click Data Review.

-Or-

Click **Data Review** in the navigation pane.

2. In the Data Review navigation pane, 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, see "Working with Master Methods" on page 88.

- 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**.

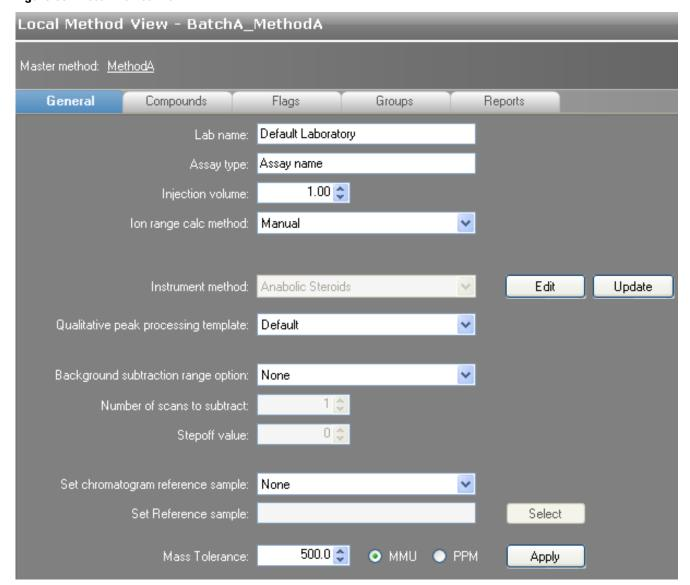


The application overwrites the local method with the master method of the same name. You can use this feature to overwrite an edited local method with the original master method or to overwrite the local method with an updated master method.

6 Using the Data Review Mode

Working in the Local Method View

Figure 85. Local Method View



Reports

This appendix contains information about standard and custom reports.

Contents

- Specifying Reports
- Report Flags
- Sample Standard Reports

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 reports listed in this appendix meet the requirements of various methods and worldwide regulatory agencies and are designed to help track the performance of the system and method. The TraceFinder application can produce both standard reports and custom reports.

Specifying Reports

As a user in the ITAdmin or LabDirector role, you can configure a list of reports that are available for the Method Development or Acquisition mode.

For detailed information about configuring reports in the Configuration mode, see "Specifying the Reports Configuration" on page 58.

For detailed information about specifying reports when you create a method in the Method Development mode, see "Editing the Reports Page" on page 173.

For detailed information about viewing batch reports in the Acquisition mode, see "Selecting and Reviewing Reports" on page 233.

A Reports Specifying Reports

Standard Reports

For each standard report you generate, you can create a version in hardcopy print, as a PDF (.pdf) file, or in an XML (.xml) output format. In addition to the report type, you can specify a report description for each of your reports. The default report description is the report name.

The TraceFinder application can generate the following types of standard reports:

- Batch Report
- Batch Report Rev 1
- Blank 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
- High Density Internal Standard Report Long
- High Density Sample Report 1
- High Density Sample Report 1 Long
- High Density Sample Report 2
- High Density Sample Report 2 Long
- High Density Sample Report 3
- High Density Sample Report 3 Long
- Internal Standard Summary Report
- Ion Ratio Failure Report
- LCSLCSD Report (Environmental and Food Safety only)
- Manual Integration Report
- Method Detection Limit Report (Environmental and Food Safety only)
- Method Report
- Method Validation Report (Environmental and Food Safety only)
- MSMSD Report (Environmental and Food Safety only)
- Quantitation Report
- Quantitation Report 2
- Solvent Blank Report
- Surrogate Recovery Report
- TIC Report
- TIC Summary Report

To view an example of each type of standard report, see "Sample Standard Reports" on page 357.

Custom Reports

For each custom report you generate, you can create a hardcopy printout or an Excel Macro-Enabled Workbook (.xlsm) output file. The default report description is the report name.

A user in the ITAdmin or LabDirector role can configure custom reports to generate a single report for an entire batch or to create separate reports for each sample. Rather than creating separate reports for each sample, this method uses data from only the last sample to create a single report for the entire batch.

The TraceFinder application includes the following custom reports:

- AltCalibrationReport
- Alternate BatchReport
- Alternate CalibrationReport
- Alternate ConfirmationReport
- Alternate MatrixSpikeReport
- Alternate SampleReport
- Alternate SummaryReport
- BatchReport
- BlankReport
- CalibrationDensityReport
- CalibrationReport
- CheckStandardReport
- CompoundCalibrationReport
- ConfirmationReport
- ConfirmationReport2
- HighDensitySampleReport1Long
- HighDensitySampleReport2Long
- HighDensitySampleReport3Long
- HighDensitySampleReport4
- HighDensitySampleReport5
- QuantitationReport
- SteroidAnalysisReport

Target Screening Reports

- Target Screening Long Report
- Target Screening Summary Report

Report Flags

When generating or viewing a report, you might see one of the following quantification or calibration flags listed on the page.

Table 89. Quantification flags

| Flag | Definition |
|--------|--|
| Ь | Compound was observed at a concentration in a Matrix Blank sample above the specified limit. |
| S | Compound was observed at a response in a solvent blank sample above the specified limit. |
| J | Compound was observed at a concentration above the limit of detection, but below the limit of quantitation. |
| I or * | Confirming/qualifying ion ratio for a compound was observed outside the target ratio range or the coelution between quantification and confirming/qualifying ion was larger than acceptable limit. |
| С | Compound was observed at a concentration above the specified carryover limit. |
| ? | Compound was observed at a concentration above the specified linearity limit. |
| D | Compound was observed at a concentration below the specified limit of detection. |
| Q | Compound was observed at a concentration below the specified limit of quantitation. |
| POS | Compound was observed at a concentration above the specified cutoff. |

Table 90. Calibration flags

| Flag | Definition |
|---------|---|
| D | Calibration for this compound exceeded the specified maximum percent relative standard deviation (%RSD). |
| F | Response factor for this compound was below the specified minimum response factor (Min RF). |
| R | Calibration for this compound was below the specified minimum correlation coefficient (r ²). |
| A | Back calculation of the calibration points for this compound exceeded the specified maximum percent difference (Max %D). |
| X | Calibration point for this compound was excluded from the overall calibration by manual selection. |
| X(ISNF) | Calibration point for this compound was excluded from the overall calibration because its associated internal standard was not found. |

A flags failure is identified by an asterisk (*), a shaded row, or the word Fail.

Values on a report that are the result of a manual integration use an uppercase M to signify a manually integrated quantification ion and a lowercase m to signify a manually integrated qualifying/confirming ion. On alternate reports, manual integration uses a black box around the value.

Sample Standard Reports

This section shows samples of the following standard report types:

- Batch Report
- Batch Report Rev 1
- Blank 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
- High Density Internal Standard Report Long
- High Density Sample Report 1
- High Density Sample Report 1 Long
- High Density Sample Report 2
- High Density Sample Report 2 Long
- High Density Sample Report 3
- High Density Sample Report 3 Long
- Internal Standard Summary Report
- Ion Ratio Failure Report
- LCSLCSD Report (Environmental and Food Safety only)
- Manual Integration Report
- Method Detection Limit Report (Environmental and Food Safety only)
- Method Report
- Method Validation Report (Environmental and Food Safety only)
- MSMSD Report (Environmental and Food Safety only)
- Quantitation Report
- Quantitation Report 2
- Solvent Blank Report
- Surrogate Recovery Report
- TIC Report
- TIC Summary Report

Tip To easily view reports in landscape format, choose **View > Rotate View > Clockwise** from the Adobe Acrobat viewer menu.

Batch Report

| | h | K | ер | or | t | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
|--------------|--------------------------|-----------------------|--------------------------|---------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|-----------------------|-----------------------|-----------------------|-----------------------|-----------------------|-----------------------|-----------------------|-----------------------|-----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|----------------------|
| | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| | | | | Comment | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| | | | | Inj Vol Conv Factor | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 | 5.000 1.0 |
| | 0441101 040644 | R1_FDF_040511 0511 | R1.calx | Vial Pos | 2 | ω | Q | Q | 2 | Q | 5 | ω | 5 | ω | 5 | ω | Q | 2 | ω | 5 | Ω | 5 | Q | 5 | 2 | 5 | 2 | 2 | co | 2 | 2 | ω | Ω | 2 | Ω | ω | Q | c) | 2 | 2 | 2 | ĸ |
| Batch Report | 1000 C | | Call File: 011HPLR1.calx | Sample Type | Matrix Blank | Cal Std | Unknown | Unknown | Unknown | Unknown | Chk Std | Unknown | Unknown | Unknown | Chk Std | Unknown | Chk Std | Unknown | Inknown |
| Batch | | | | Level | N/A | 1XLOQ | 2XLOQ | 4XLOQ | 6XLOQ | 10XLOQ | 16XLOQ | 32XLOQ | N/A | N/A | N/A | N/A | 2XLOQ | N/A | N/A | N/A | 2XLOQ | N/A | 2XLOQ | N/A | Δ/N |
| | | | | Sample Name | LLOD_HP031711 | L1XLOQ_HP031711 | L2XLOQ_HP031711 | L4XLOQ_HP031711 | L6XLOQ_HP031711 | L10XLOQ_HP031711 | L16XLOQ_HP031711 | L32XLOQ_HP031711 | МеОН | 11_0041RB | 11_0040MB | 11_0037LCMS | L2XLOQ_HP031711 | 11_0001 | 11_0002 | 11_0003 | 11_0004 | 11_0005 | 11_0006 | 11_0007 | 11_0008 | 11_0009 | 11_0010 | 11_0011 | L2XLOQ_HP031711 | 11_0012 | 11_0013 | 11_0014 | 11_0015 | 11_0016 | 11_0017 | 11_0018 | 11_0019 | 11_0020 | 11_0021 | L2XL0Q_HP031711 | 11_0022 | 11 0023 |
| | | | | Sample ID | 1 | 2 | 3 | 4 | 5 | 9 | 7 | 80 | 6 | 10 | 11 | 12 | 13 | 14 | 15 | 16 | 17 | 18 | 19 | 20 | 21 | 22 | 23 | 24 | 25 | 26 | 27 | 28 | 29 | 30 | 31 | 32 | 33 | 34 | 35 | 36 | 37 | 38 |
| | | PC | | Date/Time | 3/17/2011 5:46:30 PM | 3/17/2011 6:06:33 PM | 3/17/2011 6:26:29 PM | 3/17/2011 6:46:23 PM | 3/17/2011 7:06:17 PM | 3/17/2011 7:26:11 PM | 3/17/2011 7:46:05 PM | 3/17/2011 8:06:03 PM | 3/17/2011 8:26:01 PM | 3/17/2011 8:45:57 PM | 3/17/2011 9:05:59 PM | 3/17/2011 9:26:01 PM | 3/17/2011 9:45:57 PM | 3/17/2011 10:05:53 PM | 3/17/2011 10:25:49 PM | 3/17/2011 10:45:41 PM | 3/17/2011 11:05:35 PM | 3/17/2011 11:25:39 PM | 3/17/2011 11:45:43 PM | 3/18/2011 12:05:43 AM | 3/18/2011 12:25:37 AM | 3/18/2011 12:45:37 AM | 3/18/2011 1:05:37 AM | 3/18/2011 1:25:35 AM | 3/18/2011 1:45:29 AM | 3/18/2011 2:05:23 AM | 3/18/2011 2:25:19 AM | 3/18/2011 2:45:17 AM | 3/18/2011 3:05:13 AM | 3/18/2011 3:25:07 AM | 3/18/2011 3:45:05 AM | 3/18/2011 4:05:09 AM | 3/18/2011 4:25:11 AM | 3/18/2011 4:45:19 AM | 3/18/2011 5:05:15 AM | 3/18/2011 5:25:15 AM | 3/18/2011 5:45:11 AM | 3/18/2011 6:05:03 AM |
| de lormont | Lab Name: Inelliforation | | | File Name | LLOD_HP031711 | L1XLOQ_HP031711 | L2XL0Q_HP031711 | L4XLOQ_HP031711 | L6XLOQ_HP031711 | L10XLOQ_HP031711 | L16XLOQ_HP031711 | L32XLOQ_HP031711 | MeOH_01 | 11_0041RB | 11_0040MB | 11_0037LCMS | L2XLOQ_HP031711_01 | 11_0001 | 11_0002 | 11_0003 | 11_0004 | 11_0005 | 11_0006 | 11_0007 | 11_0008 | 11_0009 | 11_0010 | 11_0011 | L2XLOQ_HP031711_02 | 11_0012 | 11_0013 | 11_0014 | 11_0015 | 11_0016 | 11_0017 | 11_0018 | 11_0019 | 11_0020 | 11_0021 | L2XLOQ_HP031711_03 | 11_0022 | 11 0023 |

Batch Report Rev 1

| 753001 OT | 20 | | | | | 041000000000000000000000000000000000000 | | | |
|--------------------|-----------------------|-----------|------------------|--------|--------------------|---|---------------------|---------|--|
| User: TQU00637RPC | 57 37/RPC | | | - | Method: UTTHPERT_P | -K1_FDF_040511 10511 | | | |
| | L | | | J | Call File: 011HPL | 011HPLR1.calx | | | |
| File Name | Date/Time | Sample ID | Sample Name | Level | Sample Type | Vial Pos | Inj Vol Conv Factor | Comment | |
| LLOD_HP031711 | 3/17/2011 5:46:30 PM | + | LLOD_HP031711 | N/A | Matrix Blank | 2 | 5.000 1.0 | | |
| L1XLOQ_HP031711 | 3/17/2011 6:06:33 PM | 2 | L1XLOQ_HP031711 | 1XLOQ | Cal Std | 2 | 5.000 1.0 | | |
| L2XLOQ_HP031711 | 3/17/2011 6:26:29 PM | 3 | L2XLOQ_HP031711 | 2XLOQ | Cal Std | S | 5.000 1.0 | | |
| L4XLOQ_HP031711 | 3/17/2011 6:46:23 PM | 4 | L4XLOQ_HP031711 | 4XLOQ | Cal Std | 2 | 5.000 1.0 | | |
| L6XLOQ_HP031711 | 3/17/2011 7:06:17 PM | 2 | L6XLOQ_HP031711 | 6XLOQ | Cal Std | S | 5.000 1.0 | | |
| L10XLOQ_HP031711 | 3/17/2011 7:26:11 PM | 9 | L10XLOQ_HP031711 | 10XLOQ | Cal Std | S | 5.000 1.0 | | |
| L16XLOQ_HP031711 | 3/17/2011 7:46:05 PM | 7 | L16XLOQ_HP031711 | 16XLOQ | Cal Std | 2 | 5.000 1.0 | | |
| L32XLOQ_HP031711 | 3/17/2011 8:06:03 PM | 80 | L32XLOQ_HP031711 | 32XL0Q | Cal Std | 2 | 5.000 1.0 | | |
| MeOH_01 | 3/17/2011 8:26:01 PM | 6 | МеОН | N/A | Unknown | υ | 5.000 1.0 | | |
| 11_0041RB | 3/17/2011 8:45:57 PM | 10 | 11_0041RB | N/A | Unknown | 5 | 5.000 1.0 | | |
| 11_0040MB | 3/17/2011 9:05:59 PM | 11 | 11_0040MB | N/A | Unknown | υ | 5.000 1.0 | | |
| 11_0037LCMS | 3/17/2011 9:26:01 PM | 12 | 11_0037LCMS | N/A | Unknown | S | 5.000 1.0 | | |
| L2XLOQ_HP031711_01 | 3/17/2011 9:45:57 PM | 13 | L2XLOQ_HP031711 | 2XLOQ | Chk Std | S | 5.000 1.0 | | |
| 11_0001 | 3/17/2011 10:05:53 PM | 14 | 11_0001 | N/A | Unknown | c) | 5.000 1.0 | | |
| 11_0002 | 3/17/2011 10:25:49 PM | 15 | 11_0002 | N/A | Unknown | 2 | 5.000 1.0 | | |
| 11_0003 | 3/17/2011 10:45:41 PM | 16 | 11_0003 | N/A | Unknown | 2 | 5.000 1.0 | | |
| 11_0004 | 3/17/2011 11:05:35 PM | 17 | 11_0004 | N/A | Unknown | 2 | 5.000 1.0 | | |
| 11_0005 | 3/17/2011 11:25:39 PM | 18 | 11_0005 | N/A | Unknown | ις | 5.000 1.0 | | |
| 11_0006 | 3/17/2011 11:45:43 PM | 19 | 11_0006 | N/A | Unknown | 2 | 5.000 1.0 | | |
| 11_0007 | 3/18/2011 12:05:43 AM | 20 | 11_0007 | N/A | Unknown | ις | 5.000 1.0 | | |
| 11_0008 | 3/18/2011 12:25:37 AM | 21 | 11_0008 | N/A | Unknown | c) | 5.000 1.0 | | |
| 11_0009 | 3/18/2011 12:45:37 AM | 22 | 11_0009 | N/A | Unknown | 2 | 5.000 1.0 | | |
| 11_0010 | 3/18/2011 1:05:37 AM | 23 | 11_0010 | N/A | Unknown | ις | 5.000 1.0 | | |
| 11_0011 | 3/18/2011 1:25:35 AM | 24 | 11_0011 | N/A | Unknown | 2 | 5.000 1.0 | | |
| L2XLOQ_HP031711_02 | 3/18/2011 1:45:29 AM | 25 | L2XLOQ_HP031711 | 2XLOQ | Chk Std | 2 | 5.000 1.0 | | |
| 11_0012 | 3/18/2011 2:05:23 AM | 26 | 11_0012 | N/A | Unknown | 2 | 5.000 1.0 | | |
| 11_0013 | 3/18/2011 2:25:19 AM | 27 | 11_0013 | N/A | Unknown | ις | 5.000 1.0 | | |
| 11_0014 | 3/18/2011 2:45:17 AM | 28 | 11_0014 | N/A | Unknown | 2 | 5.000 1.0 | | |
| 11_0015 | 3/18/2011 3:05:13 AM | 29 | 11_0015 | N/A | Unknown | 2 | 5.000 1.0 | | |
| 11_0016 | 3/18/2011 3:25:07 AM | 30 | 11_0016 | N/A | Unknown | 2 | 5.000 1.0 | | |
| 11_0017 | 3/18/2011 3:45:05 AM | 31 | 11_0017 | N/A | Unknown | c) | 5.000 1.0 | | |
| 11_0018 | 3/18/2011 4:05:09 AM | 32 | 11_0018 | N/A | Unknown | 2 | 5.000 1.0 | | |
| 11_0019 | 3/18/2011 4:25:11 AM | 33 | 11_0019 | N/A | Unknown | 2 | 5.000 1.0 | | |
| 11_0020 | 3/18/2011 4:45:19 AM | 34 | 11_0020 | N/A | Unknown | 2 | 5.000 1.0 | | |
| 11_0021 | 3/18/2011 5:05:15 AM | 35 | 11_0021 | N/A | Unknown | 2 | 5.000 1.0 | | |
| L2XLOQ_HP031711_03 | 3/18/2011 5:25:15 AM | 36 | L2XLOQ_HP031711 | 2XLOQ | Chk Std | υ | 5.000 1.0 | | |
| 11_0022 | 3/18/2011 5:45:11 AM | 37 | 11_0022 | N/A | Unknown | S | 5.000 1.0 | | |
| 11_0023 | 3/18/2011 6:05:03 AM | 38 | 11_0023 | N/A | Unknown | 5 | 5.000 1.0 | | |

Blank Report

| b Name: | Thermo Lab | | | | | | | | | | | | Page 1 of 6 |
|-----------------|--------------|----------------------------|--------|--------------------|------------|-----------------|---------------------|---------------------|-----------------------------------|---------|----------|------|-------------|
| frument: | TQU00637 | | | | | Method: 0111 | 011HPLR1_PDP_040511 | 511 | | | | | |
| Ë | TQU00637/RPC | | | | | PDP | PDP_040511 | | | | | | |
| tch: | 011HPLR1 | | | | | Call File: 011h | 011HPLR1.calx | | | | | | |
| lal Pos | Sample ID | File Name LLOD_HP031711 | | Level N/A | | Sample Name | | File Date 3/17/2011 | File Date 3/17/2011 5:46:30 PM | Comment | Till I | | |
| | | 1 | | | | Average RF/ | Injected | 1 | Calculated | | , | | |
| Propoxur_(S) | | 7.69 | 111.10 | response 155388 | Quadratic | 000 | 5.89 | BPB 84 | 5.89 | B dd | 3.00 | 匝 | riags D |
| | | | | | | Average RF/ | petpelul | l | Calculated | l | | | |
| irget Compounds | <u>so</u> | 늏 | Olo | Response | Curve Type | Response Ratio | Conc | Units | Conc | Units | Max Conc | | Flags |
| Oxamyl_Oxime | | 4.32 | 72.10 | 1291350 | Quadratic | 0.00 | 12.67 | PPB | 12.67 | PPB | 6.00 | Tall | ٩ |
| Omethoate | | 4.34 | 183.04 | 540508 | Quadratic | 0.00 | 1.57 | PPB | 1.57 | PPB | 0.75 | 歷 | ۵ |
| Formetanate | | 4.35 | 165.10 | 4408552 | Quadratic | 0.00 | 6.27 | PPB | 6.27 | PPB | 3.00 | æ | ۵ |
| Dinotefuran | | 4.52 | 129.14 | 608336 | Quadratic | 0.00 | 6.80 | PPB | 6.80 | PPB | 3.00 | 匝 | a |
| Aldicarb_SO | | 4.51 | 132.05 | 538697 | Quadratic | 0.00 | 6:39 | PPB | 6:39 | PPB | 3.00 | Ē | ۵ |
| Propamocarb | | 4.53 | 102.00 | 750180 | Quadratic | 0.00 | 2.54 | PPB | 2.54 | PPB | 3.00 | Pass | |
| Pymetrozine | | 4.56 | 105.10 | 231305 | Quadratic | 0.00 | 1.59 | PPB | 1.59 | PPB | 0.75 | 豆 | ٩ |
| Aldicarb_SO2 | | 4.68 | 148.04 | 842061 | Quadratic | 0.00 | 6.46 | PPB | 6.46 | PPB | 3.00 | Ē | ۵ |
| Oxamyl | | 4.77 | 90.00 | 408081 | Quadratic | 0.00 | 6.44 | PPB | 6.44 | PPB | 3.00 | Ē | ۵ |
| Methomyl | | 4.98 | 88.10 | 973424 | Quadratic | 0.00 | 17.35 | PPB | 17.35 | PPB | 6.00 | Ē | a |
| Flonicamid | | 5.05 | 203.00 | 121206 | Quadratic | 0.00 | 17.91 | PPB | 17.91 | PPB | 9.00 | Ē | a |
| ODMS | | 5.07 | 169.00 | 495002 | Quadratic | 0.00 | 3.18 | PPB | 3.18 | PPB | 1.50 | Fail | a |
| 5-0H_TBZ | | 5.12 | 147.10 | 427244 | Quadratic | 0.00 | 2.75 | PPB | 2.75 | PPB | 1.50 | 屋 | ٩ |
| Thiamethoxam | | 5.12 | 181.10 | 149534 | Quadratic | 0.00 | 1.62 | PPB | 1.62 | PPB | 0.75 | Ē | ۵ |
| Monocrotophos | | 5.26 | 193.00 | 820046 | Quadratic | 0.00 | 3.22 | PPB | 3.22 | PPB | 1.50 | 匝 | a |
| Imidacloprid | | 5.66 | 209.06 | 611636 | Quadratic | 0.00 | 6.51 | PPB | 6.51 | PPB | 3.00 | Ē | a |
| Clothianidin | | 5.71 | 169.10 | 216332 | Quadratic | 0.00 | 3.14 | PPB | 3.14 | PPB | 1.50 | E E | a |
| Thiabendazole | | 5.87 | 131.10 | 692435 | Quadratic | 0.00 | 2.99 | PPB | 2.99 | PPB | 1.50 | 豆 | a |
| 3-OH_Carbofuran | an | 6.03 | 163.08 | 1717688 | Quadratic | 0.00 | 6.43 | PPB | 6.43 | PPB | 3.00 | Ē | a |

Flag legend: LOD<J<LOQ; I=lon ratio failure; C=Carryover; ?=Linearity limit; D=Detection limit; Q=Quan limit; POS=Rpt limit; b=Blank; s=Solvent blank

Calibration Report

Calibration Report

Page 1 of 7

Thermo Lab TQU00637 011HPLR1_PDP_040511 Method:

TQU00637\RPC PDP_040511 011HPLR1 011HPLR1.calx Cali File:

Calibration Summary

Lab Name:

| Calibration Summary | | | | | | | |
|---------------------|------------|------------|-------------|---------|-----------|--------|------|
| | | | A0 | A1 | A2 | R^2 | |
| | Manually | | y-Intercept | Slope | | R^2 | |
| Compound | Integrated | Curve Type | Mean RF | | | % RSD | Flag |
| Oxamyl_Oxime | | Q | 0e0 | 1.021e5 | -9.21e0 | 0.9999 | |
| Omethoate | | Q | 0e0 | 3.449e5 | -2.743e2 | 0.9999 | |
| Formetanate | | Q | 0e0 | 7.037e5 | -8.207e1 | 0.9999 | |
| Dinotefuran | | Q | 0e0 | 8.978e4 | -3.798e1 | 0.9995 | |
| Aldicarb_SO | | Q | 0e0 | 8.444e4 | -1.968e1 | 0.9997 | |
| Propamocarb | | Q | 0e0 | 2.953e5 | 1.094e2 | 0.9990 | |
| Pymetrozine | | Q | 0e0 | 1.458e5 | -1.625e2 | 0.9999 | |
| Aldicarb_SO2 | | Q | 0e0 | 1.304e5 | -2.433e1 | 0.9999 | |
| Oxamyl | | Q | 0e0 | 6.343e4 | -9.365e0 | 0.9999 | |
| Methomyl | | Q | 0e0 | 5.609e4 | 2.406e-2 | 0.9998 | A |
| Flonicamid | | Q | 0e0 | 6.78e3 | -6.227e-1 | 0.9999 | |
| ODMS | | Q | 0e0 | 1.559e5 | -3.779e1 | 1.0000 | |
| 5-OH_TBZ | | Q | 0e0 | 1.554e5 | -8.3e1 | 1.0000 | |
| Thiamethoxam | | Q | 0e0 | 9.218e4 | -8.168e1 | 0.9999 | |
| Monocrotophos | | Q | 0e0 | 2.546e5 | 3.747e0 | 0.9999 | |
| Imidacloprid | | Q | 0e0 | 9.412e4 | -1.542e1 | 0.9999 | |
| Clothianidin | | Q | 0e0 | 6.899e4 | -2.01e1 | 0.9999 | |
| Thiabendazole | | Q | 0e0 | 2.318e5 | -3.829e1 | 0.9998 | |
| 3-OH_Carbofuran | | Q | 0e0 | 2.673e5 | -5.488e1 | 0.9999 | |
| Acetamiprid | | Q | 0e0 | 1.465e5 | -1.234e2 | 0.9999 | |
| Cymoxanil | | Q | 0e0 | 1.907e4 | -2.093e0 | 0.9999 | |
| Thiacloprid | | Q | 0e0 | 1.804e5 | -3.743e1 | 0.9998 | |
| Methidathion_OA | | Q | 0e0 | 2.452e5 | -4.327e1 | 0.9999 | |
| Aldicarb | | Q | 0e0 | 9.107e4 | -8.106e0 | 0.9999 | |
| Azinphos_Me_OA | | Q | 0e0 | 7.722e4 | -1.808e1 | 0.9993 | |
| Metribuzin | | Q | 0e0 | 1.285e5 | -6.658e1 | 0.9998 | |
| Simazine | | Q | 0e0 | 5.641e4 | -4.048e1 | 0.9997 | |
| Propoxur_(S) | | Q | 0e0 | 2.644e5 | -7.519e1 | 0.9999 | |
| Pirimicarb | | Q | 0e0 | 4.722e5 | -4.082e1 | 0.9998 | |
| Bendiocarb | | Q | 0e0 | 1.491e5 | -6.818e1 | 0.9997 | |
| Carbofuran | | Q | 0e0 | 7.01e5 | -2.286e2 | 0.9998 | |
| Fenamiphos_SO | | Q | 0e0 | 1.522e4 | -9.643e-1 | 0.9998 | A |
| Tebuthiuron | | Q | 0e0 | 8.461e5 | -2.67e2 | 0.9998 | |
| Carboxin | | Q | 0e0 | 6.506e5 | -1.708e2 | 1.0000 | |
| Sulfentrazone | | Q | 0e0 | 1.374e4 | -1.836e0 | 0.9998 | |
| Fenamiphos_SO2 | | Q | 0e0 | 9.059e4 | -2.877e1 | 0.9999 | |
| Carbaryl | | Q | 0e0 | 4.251e5 | -4.87e1 | 0.9999 | |
| Thiodicarb | | Q | 0e0 | 3.968e4 | -3.693e0 | 0.9995 | A |
| Phorate_Sulfoxide | | Q | 0e0 | 3.861e5 | -1.033e2 | 0.9998 | |
| Norflurazon_DM | | Q | 0e0 | 4.987e4 | -1.219e1 | 0.9998 | |
| Phorate_Sulfone | | Q | 0e0 | 2.959e5 | -3.89e1 | 1.0000 | |
| Atrazine | | Q | 0e0 | 3.679e5 | -1.13e2 | 1.0000 | |
| Isoprocarb | | Q | 0e0 | 9.685e4 | 3.858e0 | 0.9996 | |
| Imazalil | | Q | 0e0 | 7.833e4 | -4.144e0 | 1.0000 | |
| Metalaxyl | | Q | 0e0 | 8.234e5 | 6.852e1 | 0.9999 | |
| Diuron | | Q | 0e0 | 8.495e4 | -1.208e1 | 0.9999 | |
| Norflurazon | | Q | 0e0 | 5.513e4 | -9.775e0 | 0.9999 | |
| Chlorantraniliprole | | Q | 0e0 | 2.947e4 | -1.6e0 | 0.9996 | |
| Azinphos_Me | | Q | 0e0 | 7.972e4 | -1.585e1 | 1.0000 | |
| Benoxacor | | Q | 0e0 | 1.333e4 | -8.821e-1 | 0.9999 | |
| Fluridone | | Q | 0e0 | 5.003e5 | -1.602e2 | 0.9999 | |
| Pyrimethanil | | Q | 0e0 | 4.783e4 | -1.313e0 | 0.9999 | |
| Azoxystrobin | | Q | 0e0 | 7.944e5 | -3.268e2 | 1.0000 | |

Calibration flags: D=RSD; F=Response factor; R=R^2; A=Amount;X=Excluded; X(ISNF)=Excluded because ISTD wasn't found.

A Reports

Sample Standard Reports

Calibration Report

Page 4 of 7

 Lab Name:
 Thermo Lab

 Instrument:
 TQU00637

 User:
 TQU00637\RPC

 Batch:
 011HPLR1

PDP_040511

Cali File: 011HPLR1.calx

011HPLR1_PDP_040511

| Calibration Data Points | Curve | | | | | | | |
|-------------------------|-------|----------|----------|----------|-----------|-----------|-----------|-----------|
| Compound | Туре | 1XLOQ | 2XLOQ | 4XLOQ | 6XLOQ | 10XLOQ | 16XLOQ | 32XLOQ |
| Oxamyl_Oxime | Q | 4301979 | 8239118 | 16242623 | 23873974 | 39827173 | 61070387 | 115624994 |
| Omethoate | Q | 1767322 | 3484419 | 6936118 | 10019480 | 16562654 | 25318124 | 47692071 |
| Formetanate | Q | 14839193 | 28774173 | 55703047 | 83034780 | 140328387 | 215487571 | 418192129 |
| Dinotefuran | Q | 1957462 | 3754131 | 7124741 | 10430815 | 16753892 | 24286005 | 41994838 |
| Aldicarb_SO | Q | 1800889 | 3431833 | 6655525 | 10063684 | 16448105 | 24529724 | 46055255 |
| Propamocarb | Q | 4981586 | 9714037 | 23426813 | 33421290 | 62325833 | 109908202 | 233040768 |
| Pymetrozine | Q | 827514 | 1491832 | 2850157 | 4285647 | 6910457 | 10568034 | 19185279 |
| Aldicarb_SO2 | Q | 2773066 | 5247744 | 10263769 | 15236469 | 25734192 | 39154955 | 73839209 |
| Oxamyl | Q | 1360327 | 2563147 | 4952614 | 7533418 | 12426641 | 19210270 | 36777897 |
| Methomyl | Q | 2737537 | 4939807 | 9115295 | 13452238 | 22783824 | 35397006 | 71931672 |
| Flonicamid | Q | 390564 | 802238 | 1626266 | 2324940 | 3876593 | 5915996 | 10725410 |
| ODMS | Q | 1662957 | 3110348 | 6149507 | 9205405 | 15380974 | 23845736 | 46047319 |
| 5-OH_TBZ | Q | 1563076 | 3024611 | 6173603 | 9057919 | 14758723 | 22647279 | 41244357 |
| Thiamethoxam | Q | 473839 | 910788 | 1823911 | 2662361 | 4489338 | 6790735 | 12666340 |
| Monocrotophos | Q | 2676623 | 5322359 | 10272373 | 15475446 | 25680186 | 40444172 | 81911360 |
| Imidacloprid | Q | 1961674 | 3825066 | 7339375 | 11168910 | 18451926 | 28278482 | 53957135 |
| Clothianidin | Q | 714437 | 1377702 | 2705701 | 4075866 | 6843503 | 10397614 | 20034742 |
| Thiabendazole | Q | 2395330 | 4756375 | 9238564 | 13690129 | 23391302 | 35599065 | 70334577 |
| 3-OH_Carbofuran | Q | 5631688 | 10797837 | 21032050 | 31558082 | 51838968 | 79183682 | 148721201 |
| Acetamiprid | Q | 770010 | 1489167 | 2851066 | 4289654 | 7143229 | 10810031 | 20289765 |
| Cymoxanil | Q | 377047 | 756131 | 1497988 | 2306694 | 3788308 | 5818639 | 11376041 |
| Thiacloprid | Q | 941310 | 1842377 | 3581247 | 5550042 | 9091552 | 13943056 | 27954176 |
| Methidathion_OA | Q | 5135521 | 9703625 | 18972892 | 29071643 | 47865654 | 73550478 | 139292777 |
| Aldicarb | Q | 1907088 | 3842466 | 7306696 | 11001829 | 17988401 | 28021382 | 55021609 |
| Azinphos_Me_OA | Q | 1653176 | 3209908 | 6311122 | 9387092 | 15079442 | 22160680 | 42137270 |
| Metribuzin | Q | 1382329 | 2590504 | 5034977 | 7543546 | 12429050 | 18563473 | 34333097 |
| Simazine | Q | 600586 | 1153840 | 2195023 | 3290653 | 5346202 | 7846242 | 13927839 |
| Propoxur_(S) | Q | 5511712 | 10559750 | 20690769 | 30998535 | 50521949 | 76084707 | 138563122 |
| Pirimicarb | Q | 20377224 | 38877301 | 76042738 | 114813774 | 183271544 | 281193286 | 538368936 |
| Bendiocarb | Q | 1600951 | 3142333 | 5931661 | 8910739 | 14446667 | 21691320 | 40789937 |
| Carbofuran | Q | 7314397 | 14139593 | 27712966 | 41779139 | 69378009 | 104524513 | 201180038 |
| Fenamiphos_SO | Q | 121310 | 301661 | 586312 | 869804 | 1528556 | 2435404 | 4766884 |
| Tebuthiuron | Q | 4394378 | 8642882 | 16607479 | 25289637 | 42667565 | 65032815 | 128681978 |
| Carboxin | Q | 6534810 | 12905965 | 25606866 | 38078875 | 63816950 | 99587153 | 190687947 |
| Sulfentrazone | Q | 279987 | 562967 | 1047875 | 1656952 | 2720819 | 4160971 | 8048554 |
| Fenamiphos_SO2 | Q | 837268 | 1748158 | 3488544 | 5345562 | 8905884 | 13692958 | 26046963 |
| Carbaryl | Q | 4473316 | 8624723 | 17143920 | 25672675 | 42556304 | 65943386 | 131179439 |
| Thiodicarb | Q | 490260 | 1248479 | 3213413 | 4825393 | 7837944 | 12292695 | 23881048 |
| Phorate_Sulfoxide | Q | 4159733 | 7765608 | 15382246 | 23084301 | 38335816 | 58206610 | 113125292 |
| Norflurazon_DM | Q | 1052465 | 1986723 | 3983707 | 5792912 | 9714334 | 14483066 | 26957917 |

Curve Type: A=Average RF; L=Linear; Q=Quadratic; l=Internal standard; Note: Amounts displayed for internal standards represent the ISTD Response. Calibration flags: D=RSD; F=Response factor; R=R^2; A=Amount; X=Excluded; X(ISNF)=Excluded because ISTD wasn't found.

Manually integrated

Calibration Report

 Lab Name:
 Termor Lab
 Page 7 of 7

 Instrument:
 70,00637
 Method:
 01HPLR1_PDP_040511

 User:
 70,00637
 PDP_040511

 Batch:
 011HPLR1
 Call File:
 011HPLR1.calx

| Vial Pos | Sample ID | File Name | Level | Sample Name | File Date | Comment |
|----------|-----------|------------------|--------|------------------|----------------------|---------|
| 5 | 2 | L1XLOQ_HP031711 | 1XLOQ | L1XLOQ_HP031711 | 3/17/2011 6:06:33 PM | |
| 5 | 3 | L2XLOQ_HP031711 | 2XLOQ | L2XLOQ_HP031711 | 3/17/2011 6:26:29 PM | |
| 5 | 4 | L4XLOQ_HP031711 | 4XLOQ | L4XLOQ_HP031711 | 3/17/2011 6:46:23 PM | |
| 5 | 5 | L6XLOQ_HP031711 | 6XLOQ | L6XLOQ_HP031711 | 3/17/2011 7:06:17 PM | |
| 5 | 6 | L10XLOQ_HP031711 | 10XLOQ | L10XLOQ_HP031711 | 3/17/2011 7:26:11 PM | |
| 5 | 7 | L16XLOQ_HP031711 | 16XLOQ | L16XLOQ_HP031711 | 3/17/2011 7:46:05 PM | |
| 5 | 8 | L32XLOQ_HP031711 | 32XLOQ | L32XLOQ_HP031711 | 3/17/2011 8:06:03 PM | |

Curve Type: A=Average RF; L=Linear; Q=Quadratic;l=Internal standard; Note: Amounts displayed for internal standards represent the ISTD Response.

Calibration flags: D=RSD; F=Response factor; R=R^2; A=Amount;X=Excluded; X(ISNF)=Excluded because ISTD wasn't found.

Manually integrated

Calibration Density Report

Calibration Density Report

Page 1 of 1

 Lab name:
 Thermo Fisher Laboratory

 Instrument:
 Thermo Scientific Instrument

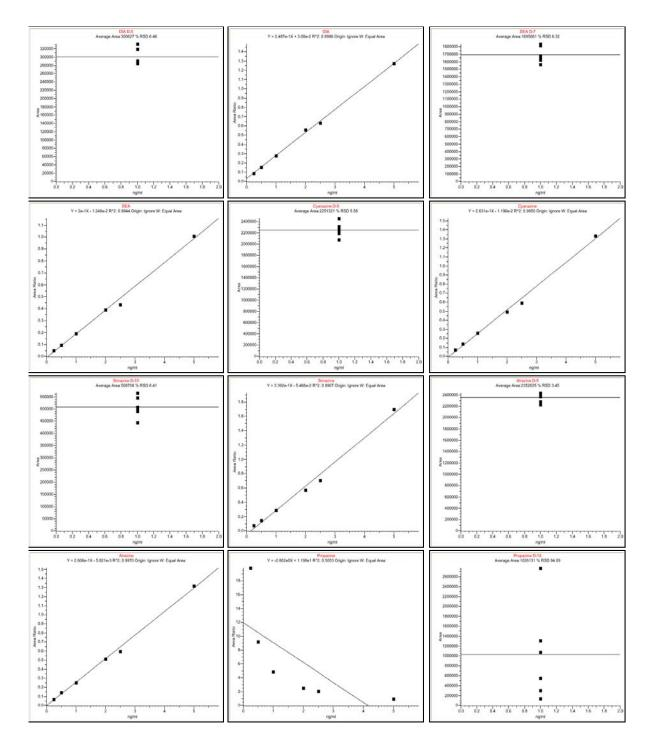
 User:
 AMER\jamie.humphries

User: AMER\jamie.humphries

Batch: Preview2

Method: Preview2_EPA536-Triazines EPA536-Triazines

Cali File: Preview2.calx



Check Standard Report

Check Standard Report

Page 1 of 4

 Lab Name:
 Thermo Lab

 Instrument:
 TQU00637

 User:
 TQU00637\RPC

011HPLR1

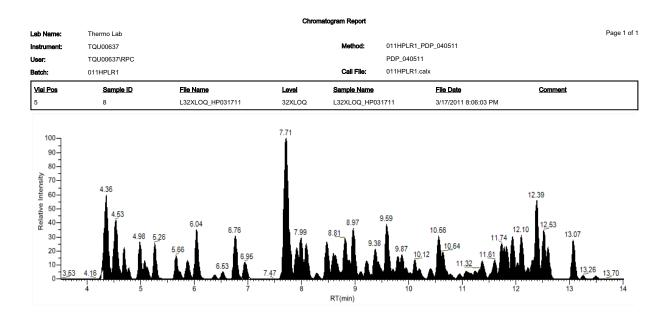
Method: 011HPLR1_PDP_040511 PDP_040511

Call File: 011HPLR1.calx

| Vial Pos | Sample ID 25 | File Name L2XLOQ_HP031711_02 | Level 2XLC | | Sample Name L2XLOQ_HP0317 | 111 | File Date 3/18/2011 1:4 | 15:29 AM | Comment | |
|-----------------|-----------------|---------------------------------|---------------|--------|---------------------------|-----------|----------------------------|----------|----------|-------------------|
| Compound | Curve Typ | | Mean RF | Min RF | RF %D | Max RF %D | QC Amt | Calc Amt | Amt %D | Max Amt %D Flag |
| Oxamyl_Oxime | Q | | | 0.000 | | | 80.00 | 79.12 | -1.11 | 20.00 Pass |
| Omethoate | Q | | | 0.000 | | | 9.91 | 9.90 | -0.08 | 20.00 Pass |
| Formetanate | Q | | | 0.000 | | | 40.10 | 39.56 | -1.34 | 20.00 Pass |
| Dinotefuran | Q | | | 0.000 | | | 40.00 | 41.32 | 3.31 | 20.00 Pass |
| Aldicarb_SO | Q | | | 0.000 | | | 40.00 | 39.04 | -2.40 | 20.00 Pass |
| Propamocarb | Q | | | 0.000 | | | 40.00 | 31.53 | -21.18 * | 20.00 Fail |
| Pymetrozine | Q | | | 0.000 | | | 10.00 | 10.38 | 3.81 | 20.00 Pass |
| Aldicarb_SO2 | Q | | | 0.000 | | | 40.20 | 39.44 | -1.88 | 20.00 Pass |
| Oxamyl | Q | | | 0.000 | | | 40.00 | 39.28 | -1.80 | 20.00 Pass |
| Methomyl | Q | | | 0.000 | | | 80.00 | 85.06 | 6.33 | 20.00 Pass |
| Flonicamid | Q | | | 0.000 | | | 120.00 | 112.98 | -5.85 | 20.00 Pass |
| ODMS | Q | | | 0.000 | | | 20.00 | 19.88 | -0.58 | 20.00 Pass |
| 5-OH_TBZ | Q | | | 0.000 | | | 20.00 | 19.03 | -4.83 | 20.00 Pass |
| Thiamethoxam | Q | | | 0.000 | | | 10.00 | 9.86 | -1.41 | 20.00 Pass |
| Monocrotophos | Q | | | 0.000 | | | 20.00 | 19.81 | -0.95 | 20.00 Pass |
| Imidacloprid | Q | | | 0.000 | | | 40.00 | 39.44 | -1.41 | 20.00 Pass |
| Clothianidin | Q | | | 0.000 | | | 20.00 | 19.84 | -0.81 | 20.00 Pass |
| Thiabendazole | Q | | | 0.000 | | | 20.00 | 19.59 | -2.04 | 20.00 Pass |
| 3-OH_Carbofuran | Q | | | 0.000 | | | 40.00 | 38.93 | -2.68 | 20.00 Pass |
| Acetamiprid | Q | | | 0.000 | | | 10.00 | 9.75 | -2.53 | 20.00 Pass |
| Cymoxanil | Q | | | 0.000 | | | 40.00 | 39.53 | -1.18 | 20.00 Pass |
| Thiacloprid | Q | | | 0.000 | | | 10.00 | 9.90 | -1.01 | 20.00 Pass |
| Methidathion_OA | Q | | | 0.000 | | | 40.00 | 39.14 | -2.15 | 20.00 Pass |
| Aldicarb | Q | | | 0.000 | | | 40.00 | 38.68 | -3.31 | 20.00 Pass |
| Azinphos_Me_OA | Q | | | 0.000 | | | 40.00 | 39.52 | -1.20 | 20.00 Pass |
| Metribuzin | Q | | | 0.000 | | | 20.00 | 19.88 | -0.59 | 20.00 Pass |
| Simazine | Q | | | 0.000 | | | 20.00 | 20.38 | 1.88 | 20.00 Pass |
| Propoxur_(S) | Q | | | 0.000 | | | 40.00 | 40.08 | 0.19 | 20.00 Pass |
| Pirimicarb | Q | | | 0.000 | | | 80.00 | 79.60 | -0.50 | 20.00 Pass |
| Bendiocarb | Q | | | 0.000 | | | 20.00 | 19.71 | -1.43 | 20.00 Pass |
| Carbofuran | Q | | | 0.000 | | | 20.00 | 19.76 | -1.19 | 20.00 Pass |
| Fenamiphos_SO | Q | | | 0.000 | | | 20.00 | 18.77 | -6.16 | 20.00 Pass |

Manually Integrated *= Fall; Curve Type: A=Average RF; L=Linear; Q=Quadratic; R=Recovery limits exceeded

Chromatogram Report



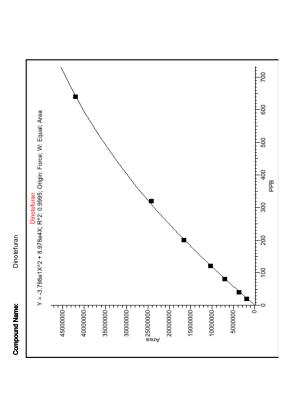
Compound Calibration Report

Page 1 of 2

011HPLR1_PDP_040511 PDP_040511 011HPLR1.calx

Method: Call File

Compound Calibration Report



| evel. | Std Amount | Std Area | IS Amount | IS Area | Response factor | Calc Amt | Units | % C/ | • |
|--------|------------|----------|-----------|---------|-----------------|----------|-------|--------|---|
| IXLOQ | 20.00 | 1957462 | | | 97873.106 | 22.01 | PPB | N/A | |
| XLOQ | 40.00 | 3754131 | | | 93853.284 | 42.58 | PPB | A/N | |
| 1XLOQ | 80.00 | 7124741 | | | 89059.260 | 82.22 | PPB | ₹/Z | |
| XLOQ | 120.00 | 10430815 | | | 86923.457 | 122.54 | PPB | ₹/Z | |
| 10XLOQ | 200.00 | 16753892 | | | 83769.461 | 204.27 | PPB | ₹/Z | |
| 16XLOQ | 320.00 | 24286005 | | | 75893.767 | 311.59 | PPB | ₹ Z | |
| 32XLOQ | 640.00 | 41994838 | | | 65616.934 | 642.34 | PPB | N/A | |
| | | | | | | | | | |
| | | | | | | | | | |
| | | | | | | | | | |

₫ ₫ ₫ ₫ ₫ Ž Ž Ž Ž Ž Ž

Thermo Scientific

Thermo Lab
TQU00637
TQU00637/RPC
011HPLR1

TraceFinder User Guide

Calibration Flags: D =RSD, F = Response factor, R = R Squared, A = Amount

| | | Compound | Compound Calibration Report | | |
|--------------|------------------|----------|-----------------------------|----------------------|-------------|
| Thermo Lab | | | | | Page 2 of 2 |
| TQU00637 | | | Method: 011HPLR1_PDP_040511 | | |
| TQU00637/RPC | | | PDP_040511 | | |
| 011HPLR1 | | | Call File 011HPLR1.calx | | |
| | | | | | |
| Sample ID | File Name | Level | Sample Name | File Date Comment | |
| 2 | L1XLOQ_HP031711 | 1XLOQ | L1XLOQ_HP031711 | 3/17/2011 6:06:33 PM | |
| 3 | L2XLOQ_HP031711 | 2XLOQ | L2XLOQ_HP031711 | 3/17/2011 6:26:29 PM | |
| 4 | L4XLOQ_HP031711 | 4XLOQ | L4XLOQ_HP031711 | 3/17/2011 6:46:23 PM | |
| 5 | L6XLOQ_HP031711 | 6XLOQ | L6XLOQ_HP031711 | 3/17/2011 7:06:17 PM | |
| 9 | L10XLOQ_HP031711 | 10XLOQ | L10XLOQ_HP031711 | 3/17/2011 7:26:11 PM | |
| 7 | L16XLOQ_HP031711 | 16XLOQ | L16XLOQ_HP031711 | 3/17/2011 7:46:05 PM | |
| 8 | L32XLOQ_HP031711 | 32XLOQ | L32XLOQ_HP031711 | 3/17/2011 8:06:03 PM | |

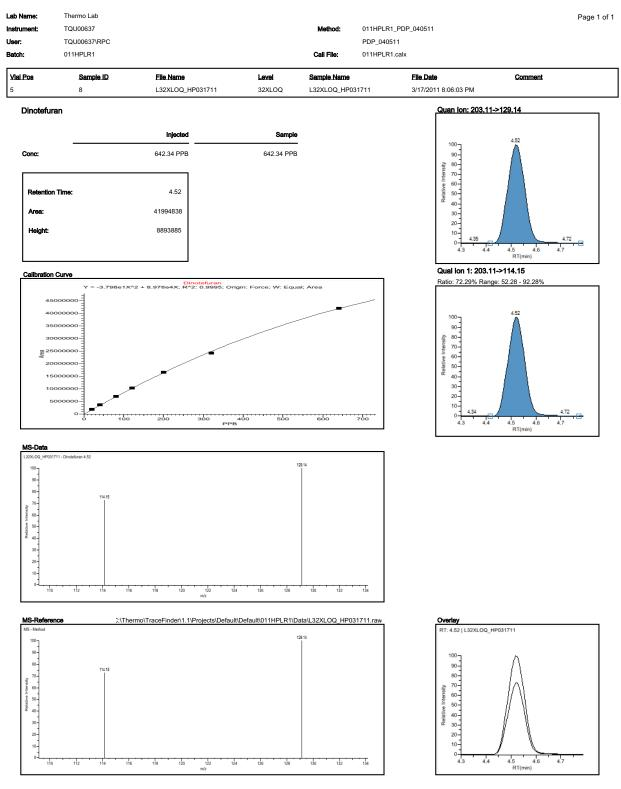
Callbration Flags: D =RSD, F = Response factor, R = R Squared, A = Amount Manually Integrated:

Compound Calibration Report - Alternate

| | | | Compound Ce | Compound Calibration Report - Alternate | | | | | Page 1 of 2 |
|---|--|----------------------------|---|--|----------|---------------------------------------|-----------|-----------|-------------|
| Lab Name: Instrument: User: Batch: | Themo Lab TQU00637 TQU00337RPC 011HPLR1 | | | Method: 011HPLR1_PDP_040511 PDP_040511 Call File 011HPLR1.calx | 040511 | | | | - NO - PBB |
| | | Compound Name: | ne: Dinotefuran | | | | | | |
| | | | Dinotefuran Y = -3.799e1X^2 + 8.978e4X; R^2: 0.9995; Origin: Force; W: Equal; Area | uran 1995; Origin: Force; W: Equal; Area | | | | | |
| | | 4500000 | 11 1111 | , | | | | | |
| | | 40000000 | | | | | | | |
| | | 35000000 | | | | | | | |
| | | 3000000 | | \ | | | | | |
| | | Area 25000000 | | \ _ | | | | | |
| | | 20000000 | | | | | | | |
| | | 1500000 | ···_ | | | | | | |
| | | 10000000 | ·•_ | | | | | | |
| | | 2000000 | • • • • • • • • • • • • • • • • • • • | | | | | | |
| | | | 0 100 200 300 | 400 500 600 | 700 | | | | |
| Quadratic | | | | | | | | | |
| Pass | File Name | Stri Area | S Amount | rea Response factor | Salo Amt | a a a a a a a a a a a a a a a a a a a | Min Range | Max Range | % RSD Flags |
| 1XLOQ | IP031711 | | | | 22.01 | В | 16.00 | 24.00 | ∀ /N |
| 2XLOQ | L2XLOQ_HP031711 40.00 | 3754131 | | 93853.284 | 42.58 | ЬРВ | 32.00 | 48.00 | N/A |
| 4XLOQ | L4XLOQ_HP031711 80.00 | 7124741 | | 89059.260 | 82.22 | PPB | 64.00 | 96.00 | Ϋ́Z |
| 6XLOQ | L6XLOQ_HP031711 120.00 | 10430815 | | 86923.457 | 122.54 | ВЬВ | 96.00 | 144.00 | N/A |
| 10XLOQ | L10XLOQ_HP031711 200.00 | 16753892 | | 83769.461 | 204.27 | ВЬВ | 160.00 | 240.00 | N/A |
| 16XLOQ | L16XLOQ_HP031711 320.00 | 24286005 | | 75893.767 | 311.59 | В | 256.00 | 384.00 | ∀/Z |
| 32XLOQ | L32XLOQ_HP031711 640.00 | 41994838 | | 65616.934 | 642.34 | PPB | 512.00 | 768.00 | Y/N |
| | | | | | | | | | |
| Manually Integrated: | grated: | Calibration Flags: D =RSD, | Calibration Flags: D =RSD, F = Response factor, R = R Squared, A = Amount | A = Amount | | | | | |

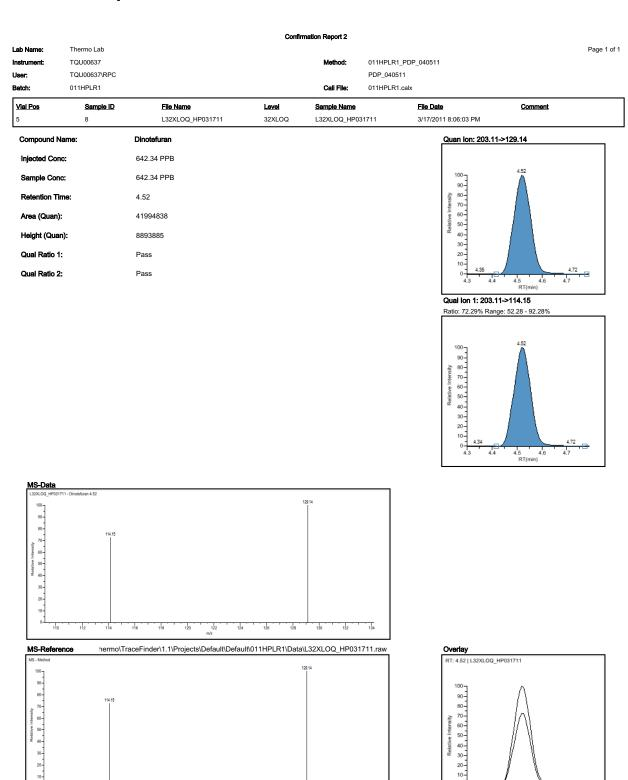
| | | | 3 | Compound Calibration Report - Attentiate | | | |
|---|--------------|------------------|--------|--|----------------------|-------------|----|
| ë | Thermo Lab | | | | | Page 2 of 2 | 01 |
| 빝 | TQU00637 | | | Method: 011HPLR1_PDP_040511 |)511 | | |
| | TQU00637/RPC | | | PDP_040511 | | | |
| | 011HPLR1 | | | Call File 011HPLR1.calx | | | |
| | | | | | | | |
| _ | Sample ID | File Name | Level | Sample Name | File Date | Comment | |
| | 2 | L1XLOQ_HP031711 | 1XL0Q | L1XLOQ_HP031711 | 3/17/2011 6:06:33 PM | | |
| | 8 | L2XLOQ_HP031711 | 2XLOQ | L2XLOQ_HP031711 | 3/17/2011 6:26:29 PM | | |
| | 4 | L4XLOQ_HP031711 | 4XLOQ | L4XLOQ_HP031711 | 3/17/2011 6:46:23 PM | | |
| | 5 | L6XLOQ_HP031711 | 6XLOQ | L6XLOQ_HP031711 | 3/17/2011 7:06:17 PM | | |
| | 9 | L10XLOQ_HP031711 | 10XLOQ | L10XLOQ_HP031711 | 3/17/2011 7:26:11 PM | | |
| | 7 | L16XLOQ_HP031711 | 16XLOQ | L16XLOQ_HP031711 | 3/17/2011 7:46:05 PM | | |
| | c | 77170000 | 000 | 771700017 | | | |

Confirmation Report

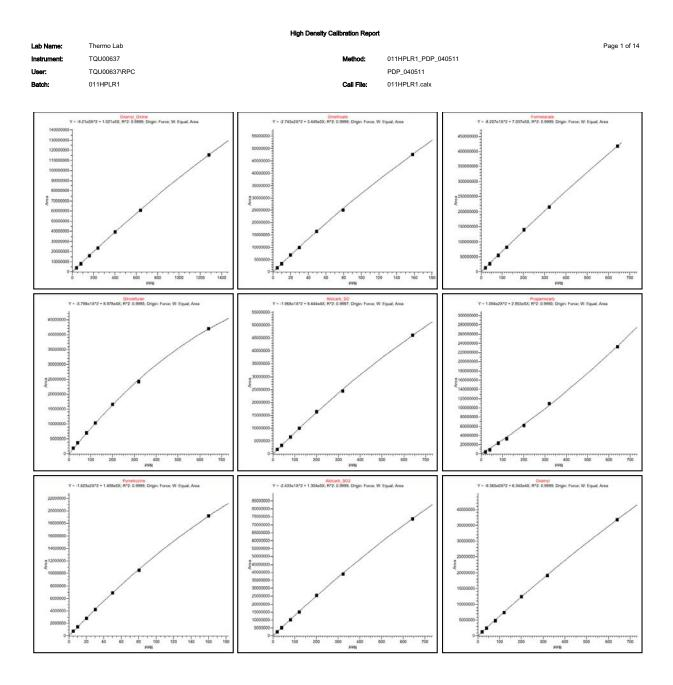


Flag legend: LOD<J<LOQ; I=Ion ratio failure; C=Carryover; ?=Linearity limit; D=Detection limit; Q=Quan limit; POS=Rpt limit; b=Blank; s=Solvent blank

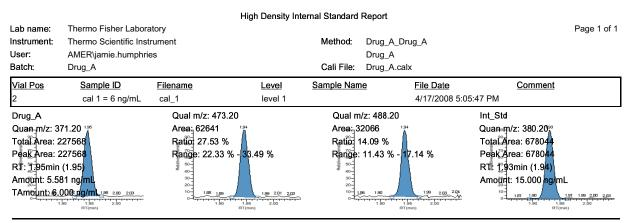
Confirmation Report 2



High Density Calibration Report



High Density Internal Standard Report



Flag legend: LOD<J<LOQ; I=Ion ratio failure; C=Carryover; ?=Linearity limit; D=Detection limit; Q=Quan limit; POS=Cutoff; n=Negative; b=Solvent blank; H=Hydrolysis

Page 1 of 1

High Density Internal Standard Report Long

High Density Internal Standard Report Long

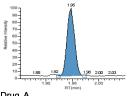
Method: Drug_A_Drug_A

Drug_A

AMER\jamie.humphries Batch: Drug_A

Cali File: Drug_A.calx

Vial Pos Sample ID <u>Filename</u> Level Sample Name File Date Comment cal 1 = 6 ng/mL cal_1 level 1 4/17/2008 5:05:47 PM



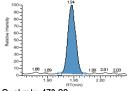
Lab name: Thermo Fisher Laboratory

User:

Instrument Thermo Scientific Instrument

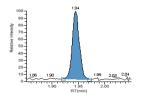
Drug_A Quan m/z: 371.20 Total Area: 227568 Peak Area: 227568

RT: 1.95min (1.95) Amount: 5.581 ng/mL TAmount: 6.000 ng/mL



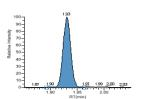
Qual m/z: 473.20 Area: 62641 Ratio: 27.53 %

Range: 22.33 % - 33.49 %



Qual m/z: 488.20 Area: 32066 Ratio: 14.09 %

Range: 11.43 % - 17.14 %

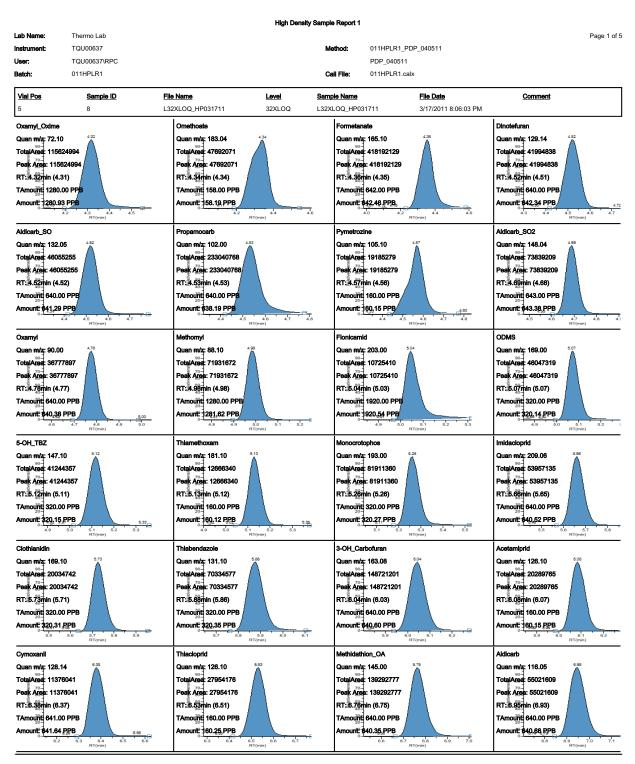


Int_Std Quan m/z: 380.20 Total Area: 678044 Peak Area: 678044 RT: 1.93min (1.94)

Amount: 15.000 ng/mL

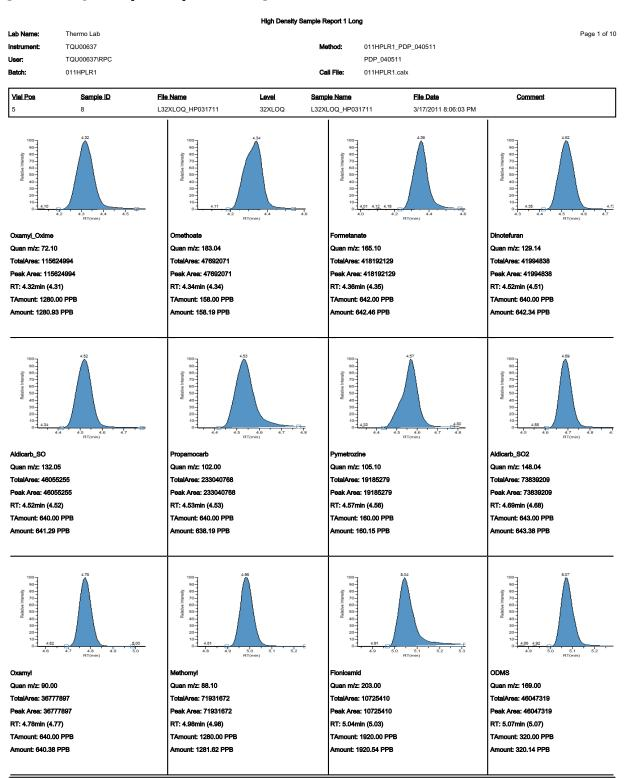
Flag legend: LOD<J<LOQ; I=lon ratio failure; C=Carryover; ?=Linearity limit; D=Detection limit; Q=Quan limit; POS=Cutoff; n=Negative; b=Solvent blank; H=Hydrolysis

High Density Sample Report 1



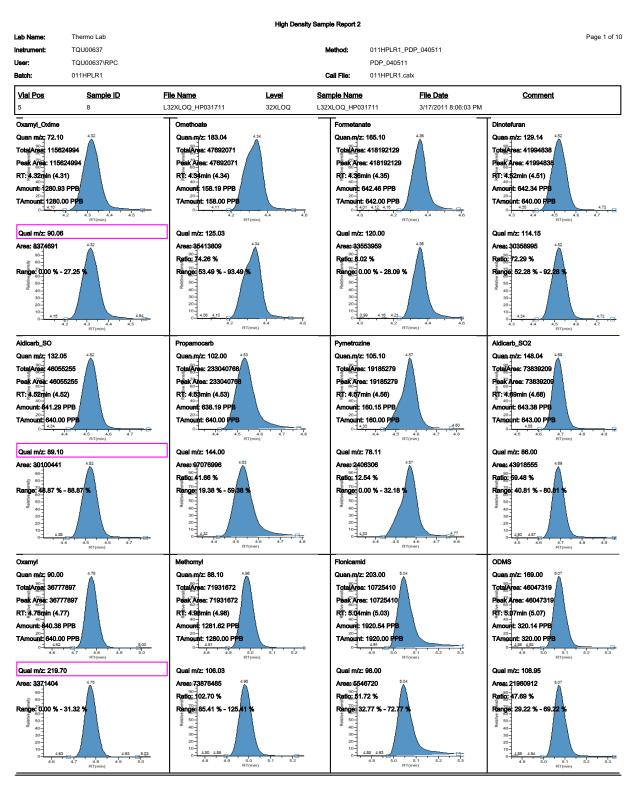
Flag legend: LOD<J<LOQ; I=Ion ratio failure; C=Carryover; ?=Linearity limit; ;D=Detection limit; Q=Quan limit; POS=Rpt limit; b=Blank; s=Solvent blank

High Density Sample Report 1 Long



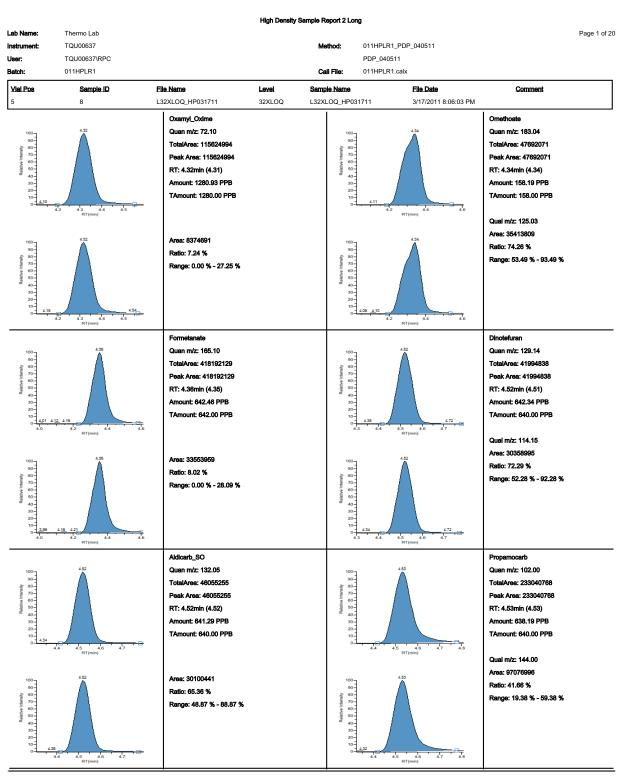
Flag legend: LOD<J<LOQ; I=Ion ratio failure; C=Carryover; ?=Linearity limit; D=Detection limit; Q=Quan limit; POS=Rpt limit; b=Blank; s=Solvent blank

High Density Sample Report 2



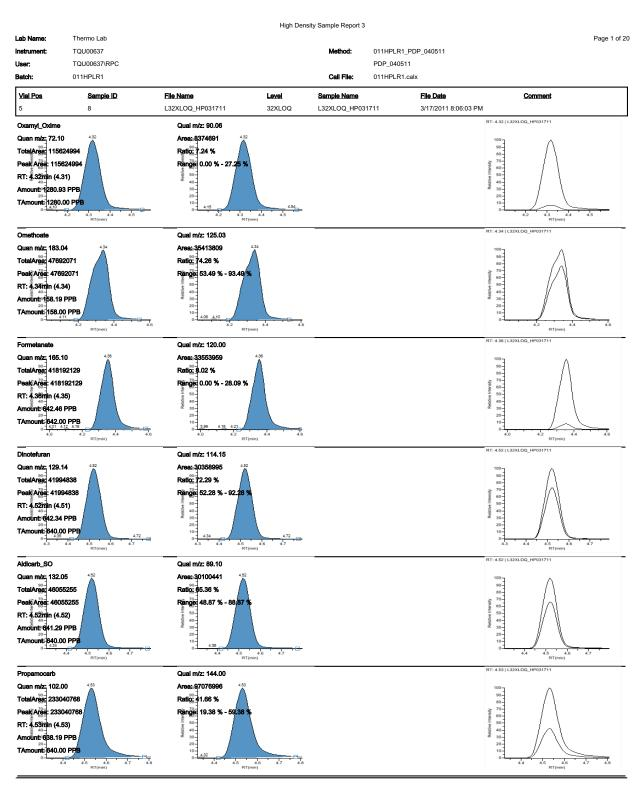
Flag legend: LOD<J<LOQ; I=Ion ratio failure; C=Carryover; ?=Linearity limit; D=Detection limit; Q=Quan limit; POS=Rot limit; b=Blank; s=Solvent blank

High Density Sample Report 2 Long



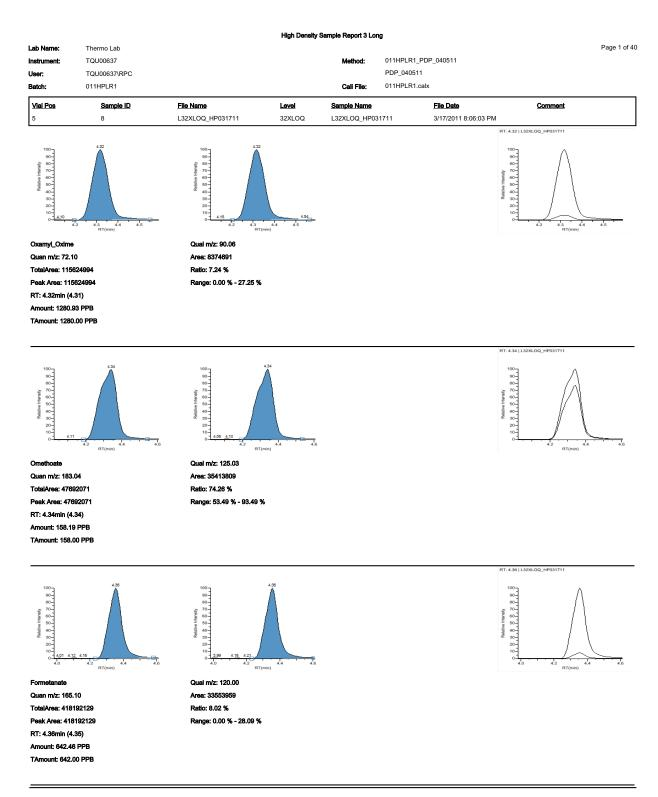
Flag legend: LOD<J<LOQ; l=lon ratio failure; C=Carryover; ?=Linearity limit; ;D=Detection limit; Q=Quan limit; POS=Rpt limit; b=Blank; s=Solvent blank

High Density Sample Report 3



Flag legend: LOD<J<LOQ; l=lon ratio failure; C=Carryover; ?=Linearity limit; D=Detection limit; Q=Quan limit; POS=Rpt limit; b=Blank; s=Solvent blank

High Density Sample Report 3 Long



Flag legend: LOD<J<LOQ; I=lon ratio failure; C=Carryover; ?=Linearity limit; D=Detection limit; Q=Quan limit; POS=Rpt limit; b=Blank; s=Solvent blank



Sample Standard Reports

Internal Standard Summary Report

Internal Standard Summary Report

Page 1 of 1

Lab name: Thermo Fisher Laboratory Method: Drug_A_Drug_A Instrument: Thermo Scientific Instrument User: AMER\jamie.humphries Drug_A Batch:

Cali File: Drug_A.calx

| Batch: | Drug_A | | | Cali File: Drug_A.o | calx | |
|-----------------|-----------------|-----------------|--------------|---------------------|----------------------|-----------------|
| <u>Vial Pos</u> | Sample ID | <u>Filename</u> | Level | Sample Name | File Date | Comment |
| 2 | cal 1 = 6 ng/mL | cal_1 | level 1 | | 4/17/2008 5:05:47 PM | |
| Compound | | | Std Response | Min | Max | Sample Response |
| Int_Std | | | 655596 | 327798(50.00%) | 983394(150.00%) | 678044 |
| | | | Std RT | Min | Max | Sample RT |
| Int Std | | | 1.93 | 1 68(-0 25) | 2 18(+0 25) | 1 93 |

* = Fail Manually integrated

Ion Ratio Failure Report

| | | | | ᅙ | Ion Ratio Fallure Report | | | | | | |
|--------------|--------------|------------------|--------|----------|--------------------------|---------------------|----------------------|---------------|--------|----------------------|---|
| ab Name: | Thermo Lab | | | | | | | | | Page 1 of 1 | _ |
| strument: | TQU00637 | | | | Method: | 011HPLR1_PDP_040511 | | | | | |
| 391 : | TQU00637/RPC | | | | | PDP_040511 | | | | | |
| atch: | 011HPLR1 | | | | Call File: | 011HPLR1.calx | | | | | |
| Vial Pos | Sample ID | File Name | Level | | Sample Name | 0 | File Date | Comment | | | _ |
| 10 | 80 | L32XLOQ_HP031711 | 32XLOQ | | L32XLOQ_HP031711 | P031711 | 3/17/2011 8:06:03 PM | | | | |
| punodwo | | | | Response | Quan lon | Quan Response | Qual lon | Qual Response | Ratio | Ratio Range | |
| sethoxydim_l | | | | Area | 178.02 | 4002074 | 220.05 | 2971396 | 74.25 | 74.25 34.09-74.09 | |
| unrofezin | | | | Δτου | 106 10 | 30861606 | 201 10 | 168167286 | 421.88 | 421 88 369 96 409 96 | |

Lab Name
Instrumen
User:
Batch:

Xial Pose
5
Compound
Sethoxydin
Buprafezin

Thermo Scientific TraceFinder User Guide

383

LCSLCSD Report (Environmental and Food Safety only)

IMPORTANT When the Sample ID is the same for an unknown sample and an LCS or LCSD sample, the unknown sample is included in the LCSLCSD report. The report information for the unknown sample displays as zeros.

| I CCI | CCD | Report |
|-------|-----|--------|
| | | |

Lab name: Thermo Fisher Laboratory Page 1 of 1

Instrument: Thermo Scientific Instrument Method: Preview2 EPA536-Triazines AMER\jamie.humphries EPA536-Triazines User:

Batch: Preview2 Cali File: Preview2.calx

| <u>Pos</u> | Sample ID | <u>Filename</u> | Level | Sample Name | File Date | Comment |
|------------|-------------|-----------------|-------|-------------|-----------------------|---------------------------|
| Tray1:16 | SampleID002 | 5ppb-002 | N/A | D008 | 6/27/2007 1:25:44 AM | New Dilutions 6/26/2007 I |
| Tray1:9 | SampleID002 | 500ppt-002 | QC | D002 | 6/26/2007 9:47:43 PM | New Dilutions 6/26/2007 I |
| Tray1:3 | SampleID002 | Cal002 | c2 | D002 | 6/26/2007 6:09:47 PM | New Dilutions 6/26/2007 I |
| Tray1:15 | SampleID008 | 5ppb-001 | N/A | D008 | 6/27/2007 12:54:35 AM | New Dilutions 6/26/2007 I |
| Tray1:8 | SampleID008 | 500ppt-001 | QC | D008 | 6/26/2007 9:16:35 PM | New Dilutions 6/26/2007 I |

| 114) 110 | Bumprens | scoppt cor | 40 | | ,000 | | 0,20,20 | 0 / 7.10.55 | | 11011 2110 | itronio o | 20,200, | |
|-------------------------|----------|------------|----------|-------|------------------|------------------|--------------|-------------|-------|------------|--------------|--------------|--|
| SampleID002 Compound | | Spike Amt | | | Lower Limit % | Upper Limit % | LCSD Conc | % Rec | RPD | Max RPD | Rec Fails | RPD Fails | |
| DIA | | 0.500 | | | 50.00 | 150.00 | 4.712 | 0.00 | 50.00 | 50.00 | 0 | 0 | |
| DEA | | 0.500 | | | 50.00 | 150.00 | 5.065 | 0.00 | 50.00 | 50.00 | 0 | 0 | |
| Cyanazine | | 0.500 | | | 50.00 | 150.00 | 5.127 | 0.00 | 50.00 | 50.00 | 0 | 0 | |
| Simazine | | 0.500 | | | 50.00 | 150.00 | 4.862 | 0.00 | 50.00 | 50.00 | 0 | 0 | |
| Atrazine | | 0.500 | | | 50.00 | 150.00 | 5.184 | 0.00 | 50.00 | 50.00 | 0 | 0 | |
| Propazine | | 0.500 | | | 50.00 | 150.00 | 3.829 | 0.00 | 50.00 | 50.00 | 0 | 0 | |
| SampleID008 | | | | | | | | | | | | | |
| Compound | | Spike Amt | LCS Conc | % Rec | Lower Limit % | Upper Limit % | LCSD Conc | % Rec | RPD | Max RPD | Rec Fails | RPD Fails | |
| DIA | | 0.500 | 4.754 | 0.00 | 50.00 | 150.00 | 4.712 | 0.00 | 0.00 | 50.00 | 0 | 0 | |
| DEA | | 0.500 | 4.960 | 0.00 | 50.00 | 150.00 | 5.065 | 0.00 | 0.00 | 50.00 | 0 | 0 | |
| Cyanazine | | 0.500 | 5.218 | 0.00 | 50.00 | 150.00 | 5.127 | 0.00 | 0.00 | 50.00 | 0 | 0 | |
| Simazine | | 0.500 | 4.839 | 0.00 | 50.00 | 150.00 | 4.862 | 0.00 | 0.00 | 50.00 | 0 | 0 | |
| Atrazine | | 0.500 | 5.178 | 0.00 | 50.00 | 150.00 | 5.184 | 0.00 | 0.00 | 50.00 | 0 | 0 | |
| Propazine | | 0.500 | 3.829 | 0.00 | 50.00 | 150.00 | 3.829 | 0.00 | 0.00 | 50.00 | 0 | 0 | |
| | | | | | | | | | | | | | |

Thermo Scientific

| Manually | integrated |
|----------|------------|
|----------|------------|

TraceFinder User Guide

384

Manual Integration Report

Manual Integration Report

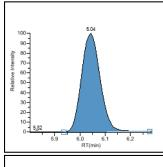
Page 1 of 2

| Lab Name: | Thermo Lab |
|-------------|--------------|
| Instrument: | TQU00637 |
| User: | TQU00637\RPC |
| Batch: | 011HPLR1 |

Method: 011HPLR1_PDP_040511
PDP_040511
Call File: 011HPLR1.calx

| | <u>fal Pos</u> | Sample ID | File Name | Level | Sample Name | File Date | Comment |
|---|----------------|-----------|------------------|--------|------------------|----------------------|---------|
| 5 | | 8 | L32XLOQ_HP031711 | 32XLOQ | L32XLOQ_HP031711 | 3/17/2011 8:06:03 PM | |

3-OH_Carbofuran m/z: 163.08

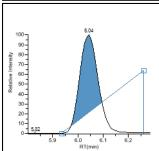


Method Integration

 Apex RT:
 6.04

 Height:
 30046630

 Area:
 148721201



Manual Integration

 Apex RT:
 6.04

 Height:
 23956657

 Area:
 94907912



Method Detection Limit Report (Environmental and Food Safety only)

Method Detection Limit Report

| Lab name: Instrument: User: Batch: | Thermo Fisher Laboratory Thermo Scientific Instrument AMER\jamie.humphries Preview2 | | | EPA | iew2_EPA536-7 536-Triazines iew2.calx | Friazines | Page 1 of 3 |
|---|---|-----------|---------|--------|---|-----------|-------------|
| Method Det | tection Limit Summary | | | | | | |
| Compound | | Avg Conc | Std Dev | t-stat | % RSD | MDL | |
| DIA D-5 | | 290218 | 0 | | 0.00 | | IS |
| DIA | | 0.095 | 0.000 | 0.000 | 0.00 | 0.000 | <<< |
| DEA D-7 | | 1704578 | 0 | | 0.00 | | IS |
| DEA | | 0.065 | 0.000 | 0.000 | 0.00 | 0.000 | <<< |
| Cyanazine | e D-5 | 2204710 | 0 | | 0.00 | | IS |
| Cyanazine | e | 0.062 | 0.000 | 0.000 | 0.00 | 0.000 | <<< |
| Simazine | D-10 | 513521 | 0 | | 0.00 | | IS |
| Simazine | | 0.168 | 0.000 | 0.000 | 0.00 | 0.000 | <<< |
| Atrazine I | D-5 | 2292164 | 0 | | 0.00 | | IS |
| Atrazine | | 0.023 | 0.000 | 0.000 | 0.00 | 0.000 | <<< |
| Propazine | | -1069.216 | 0.000 | 0.000 | 0.00 | 0.000 | <<< |
| Propazine | D-14 | 826 | 0 | | 0.00 | | IS |

Thermo Scientific

| Manually integrated | |
|---------------------|--|
| | |

TraceFinder User Guide

386

Page 2 of 3

Method Detection Limit Report

Lab name: Thermo Fisher Laboratory Instrument: Thermo Scientific Instrument $AMER \\ | jamie.humphries$ User: Batch: Preview2

Method Validation Report Data

| • | |
|----------------|-----------|
| Compound | 1 |
| DIA D-5 | 290218 |
| DIA | 0.095 |
| DEA D-7 | 1704578 |
| DEA | 0.065 |
| Cyanazine D-5 | 2204710 |
| Cyanazine | 0.062 |
| Simazine D-10 | 513521 |
| Simazine | 0.168 |
| Atrazine D-5 | 2292164 |
| Atrazine | 0.023 |
| Propazine | -1069.216 |
| Propazine D-14 | 826 |
| | |

Method: Preview2_EPA536-Triazines

EPA536-Triazines

Cali File: Preview2.calx

Manually integrated

A Reports

Sample Standard Reports

Method Detection Limit Report

Lab name: Thermo Fisher Laboratory

Method: Preview2_EPA536-Triazines

Instrument:Thermo Scientific InstrumentUser:AMER\jamie.humphries

EPA536-Triazines

Page 3 of 3

Batch: Preview2

Cali File: Preview2.calx

| Pos | Sample ID | <u>Filename</u> | Level | Sample Name | File Date | Comment |
|---------|-------------|-----------------|-------|-------------|----------------------|---------------------------|
| Tray1:1 | SampleID007 | Cal007 | N/A | D007 | 6/26/2007 8:45:28 PM | New Dilutions 6/26/2007 I |

TraceFinder User Guide Thermo Scientific

| Manually integrated | |
|---------------------|--|
|---------------------|--|

388

Page 1 of 11

Method Report

Method Report

Drug_A_Drug_A Method name:

Page number:

Master method name: Drug_A Current calibration file: Drug_A.calx

Drug_A Ion range calc method: Assay type: Average

1.000 Inj vol: Instrument method: 20x

Compound Identification:

Window View width Use as reference Reference compound Compound RT Quan mass

Int Std 380.20 1.94 10.00 0.40 Yes

0.40 No 371.20 10.00 1.95 Drug_A Int_Std

Method Report

Method name: Drug_A_Drug_A Page 2 of 11 Page number:

Master method name: Drug_A Current calibration file: Drug_A.calx

Drug_A Assay type: Ion range calc method: Average

1.000 Inj vol: Instrument method; 20x

Compound Calibration:

Compound Response Calibration Curve type Weighting Origin Units ISTD Name ISTD Units Drug_A Area Internal Linear ng/mL Int_Std ng/mL Egual Ignore

Method Report

Method name: Drug_A_Drug_A Page number: Page 3 of 11

Master method name: Drug_A Current calibration file: Drug_A.calx

Assay type: Drug_A Ion range calc method: Average

Inj vol: 1.000 Instrument method: 20x

QAQC Limits:

LÓQ Compound LÓD Cutoff ULÓL Carryover Drug_A 1.500 1.500 15.000 1000.000 1000.000

Method Report

Page 4 of 11 Drug_A_Drug_A Method name: Page number:

Master method name: Drug_A Current calibration file: Drug_A.calx

Assay type: Drug_A Ion range calc method: Average

1.000 Inj vol:

Instrument method: 20x

Groups

A Reports

Sample Standard Reports

Method Report

Average

Method name: Drug_A_Drug_A

Master method name: Drug_A

Current calibration file: Drug_A.calx

Assay type: Drug_A

Inj vol: 1.000 Instrument method: 20x

Page number:

Page 5 of 11

Report options:

Quan report options ToxLab Forms settings

Report concentration: Always Quan flags

Decimal places to be reported: Flag values below LOD: 3 True Show chromatogram on Quantitation True Flag values below LOQ: True report: Flag values above Cutoff: True Display valid compounds only False Flag values above ULOL: True

Flag values above Carryover: True True

Qual options Flag values between LOD and LOQ: Sort Qual results by: Search Index

Ion range calc method:

Enable limiting peaks: False

Calculated amount option Limit Peaks to: Calculate concentration as: Truncated

User interface options Tune time tracking options

Enable tune time tracking: False Shade row when sample is outside of False Tune file lifetime (hrs): N/A evaluation criteria:

Separate ion overlay display: True Use alternative calibration report format: False

Display quan flags and legend True

Method Report

Drug_A_Drug_A Method name: Page number: Page 6 of 11

Master method name: Drug_A Current calibration file: Drug_A.calx

Assay type: Drug_A Ion range calc method: Average

Inj vol: 1.000 Instrument method: 20x

QAQC Calibration

Compound Max RSD (%) Min RF R^2 threshold Max amt diff (%) Drug_A 20.00 0.00 0.990 20.000

Method Report

Method name: Drug_A_Drug_A Page number: Page 7 of 11

Master method name: Drug_A Current calibration file: Drug_A.calx

Assay type: Drug_A Ion range calc method: Average

Inj vol: 1.000 Instrument method: 20x

QAQC QC Check:

Compound Max RF diff (%) Min RF Drug_A 20.00 0.000

Method Report

Method name: Drug_A_Drug_A Master method name:

Drug_A

Page number:

Page 8 of 11

Current calibration file: Assay type:

Drug_A.calx Drug_A

1.000

20x

Ion range calc method:

Average

QAQC Negative:

Instrument method:

Compound Drug_A

Inj vol:

Criterion % of LOD Max value 1.500

Method Report

Drug_A_Drug_A Method name:

Master method name: Current calibration file: Drug_A Drug_A.calx

Assay type: Inj vol:

Instrument method:

Drug_A 1.000 20x

Ion range calc method: Average Page number:

0.25

Page number:

Page 9 of 11

QAQC ISTD:

Compound Int_Std

Min recovery (%) 50.00 Max recovery (%) 150.00

Min RT (-min)

Max RT (+min)

0.25

Method Report

Method name: Master method name:

Drug_A_Drug_A Drug_A Drug_A.calx

20x

Current calibration file: Assay type: Drug_A 1.000 Inj vol:

Ion range calc method:

Average

Page 10 of 11

QAQC Solvent Blank:

Instrument method:

Compound Int Std Drug_A

Method None

Quan Ion RT

Upper Limit %

0

Method Report

Method name:

Drug_A_Drug_A

Page number:

Page 11 of 11

Master method name: Current calibration file: Assay type:

Drug_A Drug_A.calx Drug_A

Ion range calc method:

Average

Inj vol: 1.000 Instrument method: 20x

QAQC Hydrolysis:

CompoundName Drug_A

EvaluationMethod Range

LowerLimit 15.000

UpperLimit 23.000

Batch:

Method Validation Report (Environmental and Food Safety only)

Method Validation Report

Lab name: Thermo Fisher Laboratory Page 1 of 3

 Instrument:
 Thermo Scientific Instrument
 Method:
 Preview2_EPA536-Triazines

 User:
 AMER\jamie.humphries
 EPA536-Triazines

Preview2 Cali File: Preview2.calx

Method Validation Summary

| Withou vandation Summary | | | | | | | |
|--------------------------|----------|-----------|--------|----------|----------|------|-----------|
| Compound | Avg Conc | Theo Conc | % Diff | Min Conc | Max Conc | | Max % RSD |
| DIA D-5 | 295652 | | | | | 0.00 | IS |
| DIA | 0.358 | 0.500 | -28.34 | 0.250 | 0.750 | 0.00 | 20.00 |
| DEA D-7 | 1778658 | | | | | 0.00 | IS |
| DEA | 0.602 | 0.500 | 20.45 | 0.250 | 0.750 | 0.00 | 20.00 |
| Cyanazine D-5 | 2224244 | | | | | 0.00 | IS |
| Cyanazine | 0.565 | 0.500 | 12.90 | 0.250 | 0.750 | 0.00 | 20.00 |
| Simazine D-10 | 505462 | | | | | 0.00 | IS |
| Simazine | 0.607 | 0.500 | 21.49 | 0.250 | 0.750 | 0.00 | 20.00 |
| Atrazine D-5 | 2334865 | | | | | 0.00 | IS |
| Atrazine | 0.512 | 0.500 | 2.46 | 0.250 | 0.750 | 0.00 | 20.00 |
| Propazine | 0.757 | 0.500 | 51.41 | 0.250 | 0.750 | 0.00 | 20.00 <<< |
| Propazine D-14 | 272050 | | | | | 0.00 | IS |
| | | | | | | | |

Manually integrated <= Failure

Page 2 of 3

Method Validation Report

 Lab name:
 Thermo Fisher Laboratory

 Instrument:
 Thermo Scientific Instrument

 User:
 AMER\jamie.humphries

Batch: Preview2

Method Validation Report Data

| Compound | 1 |
|----------------|---------|
| DIA D-5 | 295652 |
| DIA | 0.358 |
| DEA D-7 | 1778658 |
| DEA | 0.602 |
| Cyanazine D-5 | 2224244 |
| Cyanazine | 0.565 |
| Simazine D-10 | 505462 |
| Simazine | 0.607 |
| Atrazine D-5 | 2334865 |
| Atrazine | 0.512 |
| Propazine | 0.757 |
| Propazine D-14 | 272050 |

rection variation report

Method: Preview2_EPA536-Triazines EPA536-Triazines

Cali File: Preview2.calx

Manually integrated <= Failure



Sample Standard Reports

Method Validation Report

Lab name: Thermo Fisher Laboratory

Method: Preview2_EPA536-Triazines

Instrument: Thermo Scientific Instrument
User: AMER\jamie.humphries

EPA536-Triazines

Page 3 of 3

Batch: Preview2

Cali File: Preview2.calx

| Pos | Sample ID | <u>Filename</u> | Level | Sample Name | File Date | Comment |
|----------|-------------|-----------------|-------|-------------|-----------------------|---------------------------|
| Tray1:10 | SampleID003 | 500ppt-003 | N/A | D003 | 6/26/2007 10:18:49 PM | New Dilutions 6/26/2007 I |

Manually integrated <= Failure

MSMSD Report (Environmental and Food Safety only)

MSMSD Report

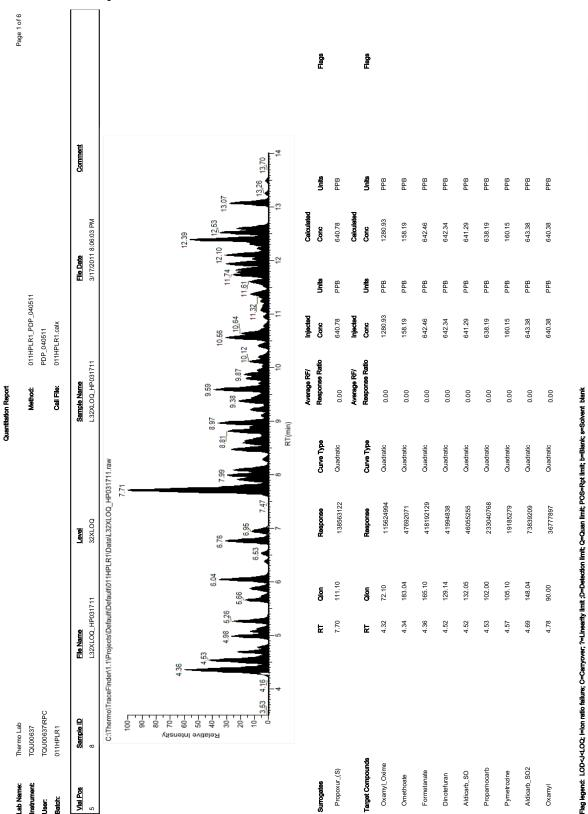
Page 1 of 1 Lab name: Thermo Fisher Laboratory Instrument: Thermo Scientific Instrument Method: Preview2_EPA536-Triazines

AMER\jamie.humphries EPA536-Triazines

User: Batch: Cali File: Preview2.calx

| Pos Tray1:1 | Sample ID SampleID021 | Filename DACTTest001 | | <u>Level</u> N/A | San D02 | nple Name | | File Date 6/27/2007 11:42:07 AM | Comment New dilution of DACT |
|----------------------|-----------------------|-------------------------|--------------|---------------------|------------|------------------|------------------|------------------------------------|------------------------------|
| SampleID021 Compound | | Unknown Conc | Spike Amt | MS Conc | % Rec | Lower Limit % | Upper Limit % | | Rec Fails |
| DIA | | 0.000 | 0.500 | 0.000 | 0.00 | 50.00 | 150.00 | | 0 |
| DEA | | 0.000 | 0.500 | 0.000 | 0.00 | 50.00 | 150.00 | | 0 |
| Cyanazine | | 0.000 | 0.500 | 0.000 | 0.00 | 50.00 | 150.00 | | 0 |
| Simazine | | 0.000 | 0.500 | 0.000 | 0.00 | 50.00 | 150.00 | | 0 |
| Atrazine | | 0.000 | 0.500 | 0.000 | 0.00 | 50.00 | 150.00 | | 0 |
| Propazine | | 0.000 | 0.500 | 0.000 | 0.00 | 50.00 | 150.00 | | 0 |

Quantitation Report



Quantitation Report - 2

| MS Integ Params: | PDP_040511 | | | | Quant Time: | | 4/19/2011 3:39:05 PM | | | |
|---|--|---------------------|----------------|------------------------|-------------------------------------|-------|---|--------------------------|--|------------|
| Quant Method: | 011HPLR1_PDP_040511 | | | | Data File: | | L32XLOQ_HP031711 | | | |
| Title: | Assay | | | | Acq On: | | 3/17/2011 8:06:03 PM | | | |
| Last Update: | 6/2/2011 9:47:56 AM | | | | Sample: | | L32XLOQ_HP031711 | | | |
| Data Acq Method: | PDP_010411 | | | | Comment: | | | | | |
| Operator: | TQU00637/RPC | | | | Vial: | | 2 | | | |
| Instr: | TQU00637 | | | | Multiplr: | | 1.000 | | | |
| Response Via: | 011HPLR1.calx | | | | Quant Results File: | v | C:\Thermo\TraceFinde OQ_HP031711.rsx | ı¹\1.1\Projects∖Default¹ | C:\Thermo\TraceFinden1.1\Projects\Default\Default\011HPLR\\Data\L32XL\\ OQ_Hp03171.rsx | |
| Surrogates | | | | | | | | | | |
| Compound Name | | RT | Qlon | Response | Conc | Units | Dev (min) | Spike Amt | Recovery | Flags |
| 28 Propoxur_(S) | 7.70 | | 111.10 | 138563122 | 640.78 PPB | | 0.01 | 40.00 | 1601.95 | |
| Target Compounds | | | | | | | | | | |
| Compound Name | | R | Olon | Response | Conc | Units | Dev (min) | | | Flags |
| 1 Oxamyl_Oxime | 4.32 | 2 | 72.10 | 115624994 | 1280.93 PPB | | 0.01 | | | |
| 2 Omethoate | 45.4 | 4 | 183.04 | 47692071 | 158.19 PPB | | 0.01 | | | |
| 3 Formetanate | 4.36 | 9 | 165.10 | 418192129 | 642.46 PPB | | 0.01 | | | |
| 4 Dinotefuran | 4.52 | 2 | 129.14 | 41994838 | 642.34 PPB | | 0.01 | | | |
| 5 Aldicarb_SO | 4.52 | 2 | 132.05 | 46055255 | 641.29 PPB | | 0.00 | | | |
| 6 Propamocarb | 4.53 | 8 | 102.00 | 233040768 | 638.19 PPB | | 00:00 | | | |
| 7 Pymetrozine | 4.57 | 7 | 105.10 | 19185279 | 160.15 PPB | | 0.01 | | | |
| 8 Aldicarb_SO2 | 4.69 | 6 | 148.04 | 73839209 | 643.38 PPB | | 0.01 | | | |
| 9 Oxamyl | 4.78 | 80 | 90.00 | 36777897 | 640.38 PPB | | 0.01 | | | |
| 10 Methomyl | 86.4 | 80 | 88.10 | 71931672 | 1281.62 PPB | | 0.01 | | | |
| 11 Flonicamid | 5.04 | | 203.00 | 10725410 | 1920.54 PPB | | 0.01 | | | |
| 12 ODMS | 5.07 | | 169.00 | 46047319 | 320.14 PPB | | 00:00 | | | |
| 13 5-OH_TBZ | 5.12 | | 147.10 | 41244357 | 320.15 PPB | | 0.01 | | | |
| 14 Thiamethoxam | 5.13 | | 181.10 | 12666340 | 160.12 PPB | | 0.01 | | | |
| 15 Monocrotophos | 5.26 | 9 | 193.00 | 81911360 | 320.27 PPB | | 0.01 | | | |
| 16 Imidacloprid | 99.6 | | 209.06 | 53957135 | 640.52 PPB | | 0.01 | | | |
| 17 Clothianidin | 5.73 | 8 | 169.10 | 20034742 | 320.31 PPB | | 0.01 | | | |
| 18 Thiabendazole | 5.88 | | 131.10 | 70334577 | 320.35 PPB | | 0.01 | | | |
| 19 3-OH_Carbofuran | 6.04 | 4 | 163.08 | 94907912 | 537.87 PPB | | 0.01 | | | - |
| | | | | | | | | | | |
| 132XI OO HP031711 | | | | Page 7 | Page 1 of 65 6/2/2011 11:40:50AM | | | | | PDP 040511 |
| Flag legend: LOD <j<loq; i="lon</th"><td>Flag legend: LOC-J<loq, ?="Linearity" b="Blank;" blank<="" c="Carryover;" d="Detection" dos="Rpt" failure;="" i="lon" limit;="" q="Quan" ratio="" s="Solvent" td=""><td>ction limit; Q=Quan</td><td>limit; POS=Rpt</td><td>limit; b=Blank; s=Solv</td><td>ent blank</td><td></td><td></td><td></td><td>Manually integrated</td><td></td></loq,></td></j<loq;> | Flag legend: LOC-J <loq, ?="Linearity" b="Blank;" blank<="" c="Carryover;" d="Detection" dos="Rpt" failure;="" i="lon" limit;="" q="Quan" ratio="" s="Solvent" td=""><td>ction limit; Q=Quan</td><td>limit; POS=Rpt</td><td>limit; b=Blank; s=Solv</td><td>ent blank</td><td></td><td></td><td></td><td>Manually integrated</td><td></td></loq,> | ction limit; Q=Quan | limit; POS=Rpt | limit; b=Blank; s=Solv | ent blank | | | | Manually integrated | |
| | | | | | | | | | , , | |

Quantitation Report - 2

| Nome | 10 | C | Doenoneo | | | |
|-----------------|-------|--------|-----------|---------------------|-------|------------|
| ייישווים | Ž | 5 | peliodeau | 200 | (min) | effet. |
| Pyrimethanil | 9.31 | 107.00 | 15176188 | 320.09 PPB | 0.02 | |
| Azoxystrobin | 9.38 | 372.10 | 118737360 | 160.00 PPB | 0.01 | |
| Fenobucarb | 9.43 | 95.00 | 68157960 | 320.10 PPB | 0.01 | |
| Linuron | 9.48 | 160.10 | 31079975 | 639.90 PPB | 0.02 | |
| Dimethomorph_I | 9.55 | 301.10 | 8049921 | 92.89 PPB | 0.00 | |
| Fenamidone | 9.58 | 92.10 | 11883329 | 320.00 PPB | 0.01 | |
| Methiocarb | 62.6 | 169.00 | 188527803 | 644.38 PPB | 0.01 | |
| Boscalid | 0.70 | 307.10 | 16765638 | 320.43 PPB | 0.01 | |
| Fludioxonil | 9.72 | 158.00 | 11536347 | 640.67 PPB | 0.02 | |
| Mandipropamid | 9.70 | 328.21 | 36436634 | 320.31 PPB | 0.00 | |
| Sethoxydim_I | 9.74 | 178.02 | 4002074 | 22.38 PPB | 0.02 | _ |
| Promecarb | 08'6 | 109.02 | 72069251 | 320.27 PPB | 0.01 | |
| Flutolanil | 9.81 | 262.00 | 59962817 | 160.05 PPB | 0.01 | |
| Fluopicolide | 9.83 | 173.00 | 29652715 | 160.16 PPB | 0.01 | |
| Methoxyfenozide | 78.6 | 149.00 | 68163623 | 319.95 PPB | 0.01 | |
| Dimethomorph_II | 88.6 | 301.10 | 24249261 | 227.43 PPB | 0.00 | |
| Clethodim_I | 9.94 | 164.09 | 30569148 | 512.02 PPB | 0.02 | |
| Triadimefon | 9.95 | 197.00 | 34892978 | 640.89 PPB | 0.01 | |
| lmiprothrin_l | 10.02 | 151.10 | 9384265 | 68.41 PPB | 0.01 | |
| Myclobutanil | 10.03 | 70.10 | 20569388 | 639.97 PPB | 0.02 | |
| Bifenazate | 10.12 | 170.10 | 47077629 | 319.44 PPB | 0.01 | |
| Triadimenol | 10.16 | 70.00 | 17384509 | 1924.45 PPB | 0.02 | |
| Spirotetramat | 10.20 | 216.15 | 28050091 | 159.99 PPB | 0.01 | |
| lmiprothrin_II | 10.22 | 123.09 | 10608234 | 248.90 PPB | 0.01 | - |
| Cyazofamid | 10.41 | 108.00 | 50568848 | 642.61 PPB | 0.03 | |
| Fenbuconazole | 10.54 | 125.00 | 26717669 | 640.05 PPB | 0.02 | |
| Chlorpyrifos_OA | 10.56 | 278.00 | 125363774 | 641.05 PPB | 0.02 | |
| Uniconazole | 10.61 | 125.08 | 9243171 | 2559.62 PPB | 0.02 | |
| Diflubenzuron | 10.62 | 158.02 | 27577198 | 1280.31 PPB | 0.03 | |
| Tebufenozide | 10.64 | 133.00 | 114645763 | 320.30 PPB | 0.03 | |
| Flubendiamide | 10.71 | 408.07 | 20125339 | 320.06 PPB | 0.02 | |
| | | | | Page 3 of 65 | | |
| P031711 | | | 6/2/2011 | AAA 40 44 4400.0000 | | 113000 AUG |
| | | | | INALC:04:11 | | 5 - |

Page 1 of 4

Solvent Blank Report

011HPLR1

Instrument:

| | Report |
|--|--------|
| | |
| | |

011HPLR1_PDP_040511 Method:

TQU00637 TQU00637\RPC PDP_040511

Cali File: 011HPLR1.calx

| Vial Pos | Sample ID | File Name | Level | Sample Name | File Date | Comment |
|----------|-----------|---------------|-------|---------------|----------------------|---------|
| 5 | 1 | LLOD_HP031711 | N/A | LLOD_HP031711 | 3/17/2011 5:46:30 PM | |

| Surrogates | RT | Qion | Response | Method | Upper Limit |
|------------------|------|--------|----------|--------|-------------|
| Propoxur_(S) | 7.69 | 111.10 | 1553885 | None | oppor a |
| | | | | | |
| Target Compounds | RT | Qion | Response | Method | Upper Limit |
| OxamyI_Oxime | 4.32 | 72.10 | 1291350 | None | |
| Omethoate | 4.34 | 183.04 | 540508 | None | |
| Formetanate | 4.35 | 165.10 | 4408552 | None | |
| Dinotefuran | 4.52 | 129.14 | 608336 | None | |
| Aldicarb_SO | 4.51 | 132.05 | 538697 | None | |
| Propamocarb | 4.53 | 102.00 | 750180 | None | |
| Pymetrozine | 4.56 | 105.10 | 231305 | None | |
| Aldicarb_SO2 | 4.68 | 148.04 | 842061 | None | |
| Oxamyl | 4.77 | 90.00 | 408081 | None | |
| Methomyl | 4.98 | 88.10 | 973424 | None | |
| Flonicamid | 5.05 | 203.00 | 121206 | None | |
| ODMS | 5.07 | 169.00 | 495002 | None | |
| 5-OH_TBZ | 5.12 | 147.10 | 427244 | None | |
| Thiamethoxam | 5.12 | 181.10 | 149534 | None | |
| Monocrotophos | 5.26 | 193.00 | 820046 | None | |
| Imidacloprid | 5.66 | 209.06 | 611636 | None | |
| Clothianidin | 5.71 | 169.10 | 216332 | None | |
| Thiabendazole | 5.87 | 131.10 | 692435 | None | |
| 3-OH_Carbofuran | 6.03 | 163.08 | 1717688 | None | |
| Acetamiprid | 6.07 | 126.10 | 239740 | None | |
| Cymoxanil | 6.38 | 128.14 | 108494 | None | |
| Thiacloprid | 6.51 | 126.10 | 299625 | None | |
| Methidathion_OA | 6.75 | 145.00 | 1537066 | None | |
| Aldicarb | 6.94 | 116.05 | 563453 | None | |
| Azinphos_Me_OA | 6.95 | 132.00 | 485885 | None | |
| Metribuzin | 7.65 | 187.16 | 437173 | None | |
| Simazine | 7.69 | 104.10 | 182755 | None | |
| Pirimicarb | 7.71 | 182.08 | 6037105 | None | |

| Manually integrated | |
|---------------------|--|
| | |



Sample Standard Reports

Surrogate Recovery Report

Surrogate Recovery Report

Page 1 of 1

 Lab Name:
 Thermor Lab

 Instrument:
 TQU00637
 Method:
 011HPLR1_PDP_040511

 User:
 TQU00637/RPC
 PDP_040511

 Batch:
 011HPLR1
 Call File:
 011HPLR1.calx

| Vial Pos | Sample ID | File Name | Level | Sample Name | File Date | Comment |
|----------|-----------|------------------|--------|------------------|----------------------|---------|
| 5 | 8 | L32XLOQ_HP031711 | 32XLOQ | L32XLOQ_HP031711 | 3/17/2011 8:06:03 PM | |

| Compound | Conc Added | Conc Recovered | % Recovered Limits | |
|--------------|------------|----------------|------------------------|---|
| Propoxur_(S) | 40.00 | 640.78 | 1601.95 83.00 - 121.00 | R |

Manually Integrated R=Recovery limits exceeded

TIC Report

TIC Report Page 1 of 1 Thermo Scientific Instrument jbe_test_5_jbe_test_q2_c1 MS110434\ITQ jbe_test_q2_c1 jbe_test_5 jbe_test_5.calx Vial Pos File Name Level File Date Sample ID Sample Name Comment spl07 3/15/2011 3:24:54 PM L135 20 ppm Height Flag 224271083 65488496 0.000 0.000 Compound CAS Number Entry Weight m-Methoxybenzoic acid, hexadecyl 111722113 mainlib 113559 565 984 C24H40O3 376 4-Methoxyphenyl octadecyl ether 562 981 C25H44O2 376 975 C19H11F3O5 1,3-Benzenediol, o-(2-furoyl)-o'-(2-t 0 mainlib 128603 6 556 376 Relative Intensity 60-40-376.00 362.00 378.00 220 240 260 280 300 m/z 320 340 360 380 4-Methoxyphenyl octadecyl ether 100 Relative Intensity 376.00 <u>377</u>.00 300 m/z 360 380 1,3-Benzenediol, o-(2-furoyl)-o'-(2-trifluoromethylbenzoyl)-80-Relative Intensity 60-40-20 376.00 220 300 m/z 380 240 260 280 320 340 360

Thermo Scientific TraceFinder User Guide 401

Flag legend: P = Library entry selected manually

TIC Summary Report

TIC Summary Report

Page 1 of 2

Lab Name: Default Laboratory Thermo Scientific Instrument Instrument:

Method:

jbe_test_5_jbe_test_q2_c1

jbe_test_q2_c1

User: MS110434\ITQ

Batch: jbe_test_5 Cali File: jbe_test_5.calx

| Vial Pos 4 | Sample ID 6 | File Name spl07 | <u>Level</u> N/A | Sample Name | File Date 3/15/2011 3:24:54 PM | Comment L135 20 ppm | |
|----------------------|-------------------------------|--------------------|---------------------|-------------|-----------------------------------|------------------------|------|
| nternal Standa | ırds | | | | | | |
| Internal Standard | | ISTD# | RT | Response | Injected Conc Units | Sample Conc Units | |
| | | | | | | | |
| Qualitatively-ide | entified Compounds | | | | Injected | Sample | |
| Compound | | Uses ISTD# | RT | Response | Conc Units | Conc Units | Flag |
| 8-Chloro-5-quinoli | inecarboxylic acid | | 2.78 | 566067 | 0.000 | 0.000 | |
| 8-Chloro-5-quinoli | inecarboxylic acid | | 2.90 | 164769 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | nexamethyl- | | 3.07 | 2136999 | 0.000 | 0.000 | |
| Cyclotetrasiloxane | e, octamethyl- | | 4.46 | 1139812 | 0.000 | 0.000 | |
| Benzaldehyde, 2,5 | 5-bis[(trimethylsilyl)oxy]- | | 5.44 | 10912 | 0.000 | 0.000 | |
| Benzoic acid, 2-[(t | trimethylsilyl)oxy]-, trimeth | ylsilyl ester | 5.80 | 87135 | 0.000 | 0.000 | |
| Silicic acid, diethy | l bis(trimethylsilyl) ester | | 7.18 | 14550 | 0.000 | 0.000 | |
| Quinoline, 2-chlor | o-6-methoxy-4-methyl- | | 7.74 | 10864 | 0.000 | 0.000 | |
| 1,1,1,3,5,5,5-Hept | tamethyltrisiloxane | | 8.43 | 10854 | 0.000 | 0.000 | |
| 1,1,1,3,5,5,5-Hept | tamethyltrisiloxane | | 9.40 | 10900 | 0.000 | 0.000 | |
| 1,1,1,3,5,5,5-Hept | tamethyltrisiloxane | | 11.02 | 200098 | 0.000 | 0.000 | |
| 1,1,1,3,5,5,5-Hept | tamethyltrisiloxane | | 11.02 | 27413 | 0.000 | 0.000 | |
| 1,1,1,3,5,5,5-Hept | tamethyltrisiloxane | | 11.08 | 69527 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | nexamethyl- | | 11.13 | 313964 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | nexamethyl- | | 11.33 | 35676 | 0.000 | 0.000 | |
| Padimate O | | | 11.44 | 843015 | 0.000 | 0.000 | |
| 1,1,1,3,5,5,5-Hept | tamethyltrisiloxane | | 11.52 | 13886 | 0.000 | 0.000 | |
| Cyclohexa-2,5-die | ene-1,4-dione, 2-methyl-5- | (4-morpholinyl)- | 11.91 | 413305 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | nexamethyl- | | 11.96 | 523374 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | nexamethyl- | | 12.07 | 206827 | 0.000 | 0.000 | |
| 5-Acetamido-4,7-c | dioxo-4,7-dihydrobenzofur | azan | 12.11 | 91980 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | nexamethyl- | | 12.13 | 43177 | 0.000 | 0.000 | |
| 5-Acetamido-4,7-c | dioxo-4,7-dihydrobenzofur | razan | 12.33 | 15408 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | nexamethyl- | | 12.48 | 15166 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | nexamethyl- | | 12.51 | 16599 | 0.000 | 0.000 | |
| 1,1,1,3,5,5,5-Hept | tamethyltrisiloxane | | 12.58 | 12284 | 0.000 | 0.000 | |
| 5-Acetamido-4,7-c | dioxo-4,7-dihydrobenzofur | razan | 12.81 | 16241 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | nexamethyl- | | 12.84 | 11652 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | nexamethyl- | | 12.86 | 12466 | 0.000 | 0.000 | |
| | tamethyltrisiloxane | | 12.90 | 57442 | 0.000 | 0.000 | |
| 1,1,1,3,5,5,5-Hept | tamethyltrisiloxane | | 12.97 | 44518 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | | | 12.98 | 15734 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | · | | 13.02 | 35927 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | · | | 13.04 | 45368 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | · | | 13.10 | 33807 | 0.000 | 0.000 | |
| Cyclotrisiloxane, h | · | | 13.12 | 67005 | 0.000 | 0.000 | |
| • | nexamethyl- | | 13.16 | 30617 | 0.000 | 0.000 | |

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, Sample Level, and Comment columns.

Use the following 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 non selected cell.



2. Right-click and choose **Copy Down** from the shortcut menu.

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 non selected cell.



3. Right-click and choose Fill Down from the shortcut menu.

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 TraceFinder 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:

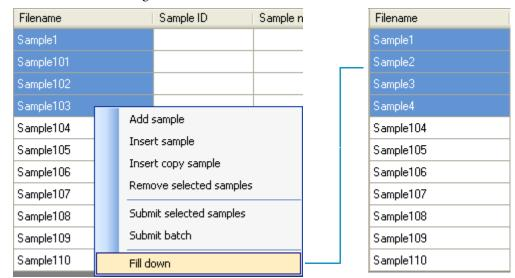
Drag your cursor to select a contiguous group of sample values.

-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.



Using Filter Criteria

The filter criteria tool is available from the compound datastore in the Configuration mode and the acquisition list in the Method Development mode.

❖ To filter the compound list

1. To display only a filtered list of compounds, click the funnel button, in the column header.

For each column, a list of filterable criteria is displayed. In all columns, your filter choices are All, Blanks, and NonBlanks. Other filter criteria are specific to the individual columns.

2. To create a custom filter based on compound values in a specific column, choose **Custom** from the column list.

The Enter Filter Criteria dialog box opens. See "Enter Filter Criteria" on page 408.

- 3. From the Operator list, select an operator.
- 4. From the Operand list, select an operand.
- 5. When all conditions are defined, click **OK**.

The complete filter string is displayed at the bottom of the dialog box, for example, chemical formula = Blanks.

Enter Filter Criteria

The Enter Filter Criteria dialog box is specifically named for the column on which you are filtering. In this example, the selected column is the Compound Name column.

Figure 86. Enter Filter Criteria dialog box

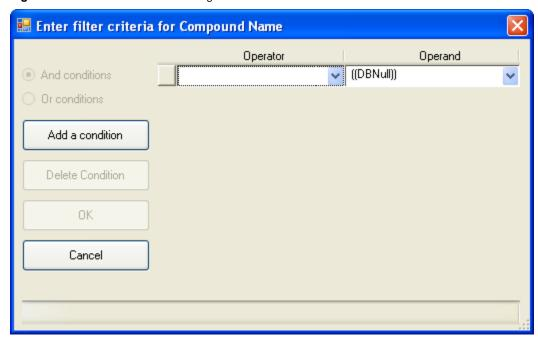


Table 91. Enter Filter Criteria dialog box parameters

| Parameter | Description |
|---------------------|---|
| And conditions | Requires meeting all filter criteria. |
| Or conditions | Requires meeting any of the specified filter criteria. |
| Add a condition | Adds a new, empty condition to the filter criteria. |
| Delete condition | Deletes the selected condition. Click the box at the left of the row to select the condition. |
| Operator | The mathematical function applied to the operand. |
| Operand | The arguments to which the operator is applied. |

Index

| Symbols | Autosampler Arm Configuration parameter /0 |
|--|--|
| csv, defined 2 | Avalon detection algorithm 71 |
| meth, defined 2 | Avalon Event List dialog box 80 |
| pmd, defined 2 | |
| raw, defined 2 | В |
| xml, defined 2 | background subtraction options 108 |
| # background scans parameter 75, 138 | Background subtraction range option parameter 112 |
| % Test parameter 155 | Baseline window parameter 77, 140 |
| 1 | Batch Level parameter 59, 176 |
| Α | Batch view 262 |
| | Batch/Template Selection view 220 |
| Account number parameter 34 | batches |
| Acquire a new raw data file parameter 96 | Acquisition mode 216 |
| Acquisition command 25 | calibration 238 |
| Acquisition List page, Master Method View 114 | Data Review mode 262 |
| Active parameter 120 | development batch 207 |
| Active View page 336 | Best match method parameter 196 |
| Add a condition parameter 408 | Blank page, Master Method View 163 |
| Add a mass as new confirming ion command 148 | Bunch Factor, event type 81 |
| Add command 80 | |
| Add compound command 52 | C |
| Add Confirming Ion command 54 | |
| Add Group command 172 | cal1-caln parameter 153 |
| Add Quan Peak command 53 | Calibration Curve page 304 |
| Add this mass to existing quan mass ranges command 148 | Calibration Levels page, Master Method View 152 |
| Add to CDS command 98 | Calibration method parameter 196 |
| Add user parameter 35 | Calibration page, Flags page, Master Method View 161 |
| Amount parameter 151 | Calibration page, Master Method View 151 |
| And conditions parameter 408 | calibration standard (Cal Std) sample type, definition 259 |
| Apply command 80 | Cancel changes parameter 35 |
| Area noise factor parameter 77, 140 | Cancel command 80 |
| Area scan window parameter 78, 141 | Carryover limit parameter 160 |
| Area tail extension parameter 78, 141 | CAS No parameter 120 |
| Area Threshold, event type 81 | Category parameter |
| Assay type parameter 111 | Acquisition List 116 Compound Datastore 48 |
| Associate a Raw Data File dialog box 99 | caution flags 289 |
| Associate a Raw Data File parameter 99 | Change command 80 |
| automated background subtraction options 108 | check standard (Chk Std) sample, definition 259 |
| Automatically create the master method parameter 96 | check standard (Chk Std) sample, definition 239 |

| Chemical Formula parameter | comma-separated values, defined 2 |
|---|--|
| Acquisition List 116 | Compound Datastore Name parameter, Save Compound |
| Compound Datastore 48 | Datastore 47 |
| Chk Std page, Master Method View 162 | Compound Datastore view 48 |
| Collision Energy parameter | Compound Datastore, administration 44 |
| Acquisition List | compound datastore, enabling 67 |
| confirming ion 118 | Compound Name parameter |
| quan peak 117 | |
| Compound Datastore | Acquisition List 116 |
| quan peak 49 | Compound Datastore 48 |
| compound datastore | Compound parameter |
| confirming ion 50 | Calibration 151 |
| commands | Calibration levels 153 |
| Acquisition 25 | Calibration page 161 |
| * | Chk Std page 162 |
| Add a mass as now son famina ion 148 | Identification 119 |
| Add a mass as new confirming ion 148 | ISTD page 166 |
| Add compound 52 | Lab Control page 168 |
| Add Confirming Ion 54 | Limits page 160 |
| Add Group 172 | Matrix Blank page 163 |
| Add Quan Peak 53 | Matrix Spike page 170 |
| Add this mass to existing quan mass ranges 148 | Method Val page 169 |
| Add to CDS 98 | QC levels page 155 |
| Apply 80 | Solvent Blank page 165 |
| Cancel 80 | Surrogate page 167 |
| Change 80 | Compound type parameter |
| Configuration 26 | Calibration 151 |
| Continue to Method 98 | Identification 120 |
| Copy down 403 | compound types |
| Data Review 25 | internal standards 119 |
| Delete 80 | quan 119 |
| Enable Target Screening 83 | surrogates 119 |
| Enable User Security 66 | target 119 |
| Exit TraceFinder 9 | Compounds page, Master Method View 114 |
| Fill down 403 | compounds, importing to compound datastore 51 |
| Help, dashboard 25 | |
| Hide Internal Standard 298 | Configuration command 26 |
| Import Compounds 51 | Configuration navigation pane 28 |
| Import Published Method 198 | Confirming Ions page 308 |
| Load Compound Datastore 46 | Confirming n Ion Ratio Flag parameter 339 |
| Log Off 25 | Confirming n Ion Ratio parameter 339 |
| Methods 26 | Confirming n Manual Flag parameter 339 |
| Modify Columns | Confirming n Mass parameter 338 |
| Batch View 269 | Confirming n Range parameter 339 |
| Data Review View 287 | Confirming n Response parameter 339 |
| New Compound Datastore 46 | Constrain peak width parameter |
| Real Time Status 25 | Genesis |
| Reset scaling 148 | |
| Save As Compound Datastore 47 | application configuration parameters 74 |
| Save Compound Datastore 46 | method development 137 |
| Set this mass as a new quan peak 148 | ICIS |
| Set this mass as quan mass 148 | application configuration 77 |
| Show Internal Standard 298 | method development 140 |
| Update confirming ion ratios with this spectrum 148 | Continue to Method command 98 |
| 1 0 | Copy down command 403 |

| Correct surrogates parameter 181 | Displayed Reports parameter |
|---|--|
| Create Master Method dialog box 89 | Environmental and Food Safety 65 |
| Create method parameter 86 | General Quantitation 62 |
| Create PDF parameter, standard reports 176 | |
| Create XLSM parameter 176 | E |
| Create XML parameter, standard reports 176 | Edit Compound Donondont Parameters dialog hov 00 |
| Curve type parameter | Edit Compound Dependent Parameters dialog box 98 |
| Calibration 151 | Edit user parameter 35 |
| Method Template Editor 196 | Email address parameter 34 |
| custom reports, list of 6 | Enable Multiplexing Support parameter 70 |
| CV Test (%) parameter 166 | Enable peak threshold parameter 196 |
| · / I | Enable Target Screening command 83 |
| D | Enable tune time tracking parameter 181 |
| ט | Enable User Security command 66 |
| dashboard parameters 25 | Enable valley detection parameter |
| Data Review command 25 | application configuration 74 |
| Data Review view 284 | method development 137 |
| Decimal places to be reported parameter 180 | Enabled parameter 34 |
| Delete command 80 | End Threshold, event type 81 |
| Delete condition parameter 408 | Energy Ramp parameter |
| Detection method parameter | application configuration 49 |
| Avalon | method development 117 |
| application configuration 79 | Event parameter 80 |
| method development 142 | Event types 81 |
| Genesis | Example parameter 176 |
| application configuration 73 | Exclude matching quan peaks parameter 198 |
| method development 137 | Exit TraceFinder command 9 |
| ICIS | Expected RT parameter, Detection 131 |
| application configuration 76 | Expected width parameter |
| method development 140 | application configuration 74 |
| Detection page, Master Method View 121 | method development 137 |
| Detector parameter, Detection 133 | Experiment Type parameter |
| Development Batch view 207 | Acquisition List 116 |
| dialog boxes | Compound Datastore 48 |
| Add Library 11 | experiment types 55 |
| Associate a Raw Data File 99 | Export SRM data parameter 87 |
| Avalon Event List 80 | Extracted Ion Chromatogram 55 |
| Create Master Method 89 | 8 |
| Edit Compound Dependent Parameters 98 | F |
| Import an Xcalibur Method 97 | Г |
| Invalid Compound Datastore not saved 46 | feature summary 3 |
| Open Chromatograph Reference Sample 109 | file types, supported 2 |
| Open Compound Datastore 46 | Fill down command 403 |
| Save Compound Datastore 47 | Filter parameter, Detection 133 |
| Select Compounds to Add 103 | Finish view 237 |
| Thermo Library Manager 11 | Flag values above Carryover parameter 181 |
| Thermo X calibur Instrument Setup 107 | Flag values above LOR parameter 181 |
| Thermo Xcalibur Roadmap 10 | Flag values above ULOL parameter 181 |
| Disable Cluster Off, event type 82 | Flag values below LOD parameter 181 |
| Disable Cluster On, event type 82 | Flag values below LOQ parameter 181 |
| Display compounds above set limit parameter 180 | Flag values between LOD and LOQ parameter 181 |
| Display Quan flags and legend parameter 180 | |

| Flags page, Master Method View 159 | Instrument method parameter |
|---|---|
| Force Cluster Off, event type 82 | General page 112 |
| Force Cluster On, event type 82 | Method Forge 96 |
| Full name parameter 34 | Instrument view 200 |
| | instruments, supported viii |
| G | Invalid Compound Datastore not saved dialog box 46 |
| | Ion Overlay page 312 |
| General page | Ion range calc method parameter 112 |
| Background subtraction range option 112 | Ionization parameter |
| editing 106 | Acquisition List 116 |
| Injection volume parameter 112 | Compound Datastore 48 |
| Instrument method parameter 112 | ISTD page, Master Method View 166 |
| Ion range calc method parameter 112 | ISTD parameter 151 |
| Number of scans to subtract parameter 112 Qualitative peak processing template parameter 112 | IT Administrator role 36 |
| Set chromatogram reference sample parameter 112 | |
| Set Reference sample parameter 112 | I |
| Stepoff value parameter 112 | L |
| General page, Master Method View 106 | Lab Control page, Master Method View 168 |
| Genesis detection algorithm 71 | lab control sample (LCS) type, defined 259 |
| | lab control sample duplicate (LCSD) type, defined 259 |
| Groups page, Master Method View 171 | Lab name parameter 111 |
| | Lens parameter |
| Н | application configuration 49 |
| Help command, dashboard 25 | method development 117 |
| Hide Internal Standard command 298 | Level parameter 155 |
| | licenses, types of ix |
| I | Limit library hits parameter 196 |
| | Limit the retention time range parameter 196 |
| ICIS detection algorithm 71 | Limits page, Master Method View 160 |
| Identification page, Master Method View 119 | Load Compound Datastore command 46 |
| Import an Xcalibur Method dialog box 97 | Local Method view 351 |
| Import Compounds command 51 | LOD (Detection limit) parameter 160 |
| Import parameter | Log Off command 25 |
| Environmental and Food Safety 65 | logging on to TraceFinder 8 |
| General Quantitation 62 | Login parameter 9 |
| Import Published Method command 198 | logon screen 9 |
| Import Xcalibur Processing Method parameter 97 | LOQ (Quantitation limit) parameter 160 |
| Include compound peak spectrum as reference spectrum | |
| parameter 196 | LOR (Reporting limit) parameter 160 |
| Include confirming ions parameter 196 | |
| Injection amount parameter 96 | M |
| Injection concentration parameter 339 | Manual Flags parameter 338 |
| Injection Units parameter 339 | Manual injection parameter 96 |
| Injection volume parameter 112 | Mass parameter |
| Installed Reports parameter | confirming ion |
| Environmental and Food Safety 65 | Acquisition List 118 |
| General Quantitation 62 | Compound Datastore 50 |
| installing | quan peak 49, 117 |
| NIST and QED libraries 10 | Master Method View |
| TraceFinder 8 | Acquisition List page 114 |
| | Blank page 163 |
| | Calibration Levels page 152 |

| Calibration page 151 | methods |
|--|--|
| Chk Std page 162 | importing Xcalibur 97 |
| Compounds page 114 | instrument 200 |
| Detection page 121 | local 351 |
| Flags page, Calibration page 161 | master 88 |
| General page 106 | Method Template Editor 195 |
| Groups page 171 | startup and shutdown 237 |
| Identification page 119 | Methods command 26 |
| ISTD page 166 | Min peak height (S/N) parameter |
| Lab Control page 168 | Genesis 75, 138 |
| Limits page 160 | ICIS 78, 141 |
| Matrix Spike page 170 | Min peak width parameter 78, 141 |
| Method Val page 169 | Min recovery (%) parameter |
| QC Levels page 154 | ISTD page 166 |
| Ratios page 149 | Lab Control page 168 |
| Real Time Viewer page 156 | Matrix Spike page 170 |
| Reports page 173 | Method Val page 169 |
| Solvent Blank page 165 | Surrogate page 167 |
| Spectrum page 144 | Min RF parameter 162 |
| Surrogate page 167 | Min RT (-min) parameter 166 |
| Master Method View, Flags page 159 | modes |
| matrix blank sample type, defined 259 | choosing 22 |
| matrix spike (MS) sample type, defined 259 | defined 25 |
| matrix spike duplicate (MSD) sample type, defined 259 | Modify Columns command |
| Matrix Spike page, Master Method View 170 | Batch View 269 |
| Max Amt Diff (%) parameter 161 | Data Review View 287 |
| Max Conc parameter 164 | Multiplet resolution parameter 78, 141 |
| Max recovery (%) parameter | multiplexing, enabling 70 |
| ISTD page 166 | maniplexing, chaomig / v |
| Lab Control page 168 | NI. |
| Matrix Spike page 170 | N |
| Method Val page 169 | Name the master method parameter 96 |
| Surrogate page 167 | Negative Peaks, event type 81 |
| Max RF Diff (%) parameter 162 | New Compound Datastore command 46 |
| Max RPD parameter | New instrument method parameter 87 |
| Lab Control page 168 | New sample list parameter 87 |
| Matrix Spike page 170 | NIST library, installing 10 |
| Max RSD (%) parameter | Noise method parameter 78, 141 |
| Calibration page 161 | Number of Scans to Subtract parameter 108 |
| Method Val page 169 | |
| Surrogate page 167 | Number of scans to subtract parameter 112 |
| Max RT (+min) parameter, ISTD page 166 | |
| method detection limits (MDL) sample type, defined 259 | 0 |
| Method parameter | Only select top peaks parameter 196 |
| Matrix Blank page 164 | Open Chromatograph Reference Sample dialog box 109 |
| Solvent Blank page 165 | Open Compound Datastore dialog box 46 |
| Method Template Editor 195 | Open instrument method parameter 87 |
| Method Val page, Master Method View 169 | Open method parameter 86 |
| | |
| method validation sample type, defined 259 | Open Qual Browser parameter 87 |
| Method view 89 | Operand parameter 408 |
| | Operator parameter 408 |
| | Or conditions parameter 408 |

| Origin parameter | Compound |
|---|---------------------------------------|
| Calibration 151 | Calibration 151 |
| Method Template Editor 197 | Calibration levels 153 |
| Overwrite parameter | Calibration page 161 |
| Method Forge dialog box 96 | Chk Std page 162 |
| Save Compound Datastore dialog box 47 | Identification 119 |
| | ISTD page 166 |
| P | Lab Control page 168 |
| r | Limits page 160 |
| parameters | Matrix Blank page 163 |
| # background scans 75, 138 | Matrix Spike page 170 |
| % Test 155 | Method Val page 169 |
| Account number 34 | QC levels page 155 |
| Acquire a new raw data file 96 | Solvent Blank page 165 |
| Active 120 | Surrogate page 167 |
| Add a condition 408 | Compound Name |
| Add user 35 | Acquisition List 116 |
| Amount 151 | Compound Datastore 48 |
| And conditions 408 | Compound type |
| Area noise factor 77, 140 | Calibration 151 |
| Area scan window 78, 141 | Identification 120 |
| Area tail extension 78, 141 | Confirming n Ion Ratio 339 |
| Assay type 111 | Confirming n Ion Ratio Flag 339 |
| Associate a Raw Data File 99 | Confirming n Manual Flag 339 |
| Automatically create the master method 96 | Confirming n Mass 338 |
| Autosampler Arm Configuration 70 | Confirming n Range 339 |
| Background subtraction range option 112 | Confirming n Response 339 |
| Baseline window 77, 140 | Constrain peak width |
| Batch Level 59, 176 | Genesis 137 |
| Best match method 196 | ICIS 77, 140 |
| cal1-caln 153 | Correct surrogates 181 |
| Calibration method 196 | Create method 86 |
| Cancel changes 35 | Create PDF, standard reports 176 |
| Carryover limit 160 | Create XLSM 176 |
| CAS No 120 | Create XML, standard reports 176 |
| Category | Curve type |
| Acquisition List 116 | Calibration 151 |
| Compound Datastore 48 | Method Template Editor 196 |
| Chemical Formula | CV Test (%) 166 |
| Acquisition List 116 | dashboard 25 |
| Compound Datastore 48 | Decimal places to be reported 180 |
| Collision Energy | Delete condition 408 |
| Acquisition List | Detection method |
| confirming ion 118 | Avalon 79, 142 |
| quan peak 117 | Genesis 73, 137 |
| compound datastore | ICIS 76, 140 |
| confirming ion 50 | Detector 133 |
| quan peak | Display compounds above set limit 180 |
| Compound Datastore 49 | Display Quan flags and legend 180 |
| | Displayed Reports |
| | Environmental and Food Safety 65 |
| | General Quantitation 62 |
| | Edit user 35 |
| | Email address 34 |

| Enable Multiplexing Support 70 | Lens |
|---|------------------------------------|
| Enable peak threshold 196 | application configuration 49 |
| Enable tune time tracking 181 | method development 117 |
| Enable valley detection | Level 155 |
| application configuration 74 | Limit library hits 196 |
| method development 137 | Limit the retention time range 196 |
| Enabled 34 | LOD (Detection limit) 160 |
| Energy Ramp | Login 9 |
| application configuration 49 | LOQ (Quantitation limit) 160 |
| method development 117 | LOR (Reporting limit) 160 |
| Event 80 | Manual Flags 338 |
| Example 176 | Manual injection 96 |
| Exclude matching quan peaks 198 | Mass |
| Expected RT, Detection 131 | confirming ion |
| Expected width | Acquisition List 118 |
| application configuration 74 | Compound Datastore 50 |
| method development 137 | quan peak 49, 117 |
| Experiment Type | Max Amt Diff (%) 161 |
| Acquisition List 116 | Max Conc 164 |
| Compound Datastore 48 | Max recovery (%) |
| Export SRM data 87 | ISTD page 166 |
| Filter, Detection 133 | Lab Control page 168 |
| Flag values above Carryover 181 | Matrix Spike page 170 |
| Flag values above LOR 181 | Method Val page 169 |
| Flag values above ULOL 181 | Surrogate page 167 |
| Flag values below LOD 181 | Max RF Diff (%) 162 |
| Flag values below LOQ 181 | Max RPD |
| Flag values between LOD and LOQ 181 | Lab Control page 168 |
| Full name 34 | Matrix Spike page 170 |
| Import | Max RSD (%) |
| Environmental and Food Safety 65 | Calibration page 161 |
| General Quantitation 62 | Method Val page 169 |
| Import Xcalibur Processing Method 97 | Surrogate page 167 |
| Include compound peak spectrum as reference | Max RT (+min), ISTD page 166 |
| spectrum 196 | Method |
| Include confirming ions 196 | Matrix Blank page 164 |
| Injection amount 96 | Solvent Blank page 165 |
| Injection concentration 339 | Min peak height (S/N) |
| Injection Units 339 | Genesis 75, 138 |
| Injection volume 112 | ICIS 78, 141 |
| Installed Reports | Min peak width 78, 141 |
| Environmental and Food Safety 65 | Min recovery (%) |
| General Quantitation 62 | ISTD page 166 |
| Instrument method | Lab Control page 168 |
| General page 112 | Matrix Spike page 170 |
| Method Forge 96 | Method Val page 169 |
| Ion range calc method 112 | Surrogate page 167 |
| Ionization | Min RF 162 |
| Acquisition List 116 | Min RT (-min) 166 |
| Compound Datastore 48 | Multiplet resolution 78, 141 |
| ISTD 151 | Name the master method 96 |
| Lab name 111 | New instrument method 87 |
| | New sample list 87 |
| | Noise method 78, 141 |

| Number of Scans to Subtract 108 | Report Title, standard reports 176 |
|--|---|
| Number of scans to subtract 112 | Response 339 |
| Only select top peaks 196 | Response via |
| Open instrument method 87 | Calibration 151 |
| Open method 86 | Method Template Editor 197 |
| Open Qual Browser 87 | Retention Time 339 |
| Operand 408 | RMS 78, 141 |
| Operator 408 | Role 34 |
| Or conditions 408 | RT |
| Origin | Acquisition List 117 |
| Calibration 151 | Calibration 151 |
| Method Template Editor 197 | Calibration levels 153 |
| Overwrite | Calibration page 161 |
| Method Forge dialog box 96 | Chk Std page 162 |
| Save Compound Datastore dialog box 47 | Compound Datastore 49 |
| Password | Identification 119 |
| logon screen 9 | ISTD page 166 |
| user administration 34 | Lab Control page 168 |
| Path 96 | Limits page 160 |
| Peak height (%) | Matrix Blank page 163 |
| Genesis 74, 137 | Matrix Spike page 170 |
| ICIS 77, 140 | Method Val page 169 |
| Peak noise factor 77, 140 | QC levels page 155 |
| Peak S/N cutoff 75, 138 | Solvent Blank page 165 |
| Percentage 164 | Surrogate page 167 |
| Phone number 34 | S/N threshold 74, 137 |
| Polarity | Sample comment 96 |
| Acquisition List 117 | Sample Concentration 339 |
| Compound Datastore 49 | Sample file 337 |
| Precursor Mass | Sample Units 339 |
| Acquisition List 116–117 | Save changes 35 |
| Compound Datastore 49–50 | Select a custom template 96 |
| Print, standard reports 176 | Select a report 337 |
| Product Mass | Select batch location 87 |
| Acquisition List 116–117 | Select Compounds from CDS 103 |
| Compound Datastore 49–50 | Sensitivity |
| QAQC Flags 338 | Avalon 79 |
| QC1-QCn 155 | Genesis 73, 137 |
| QIon 339 | ICIS 76, 140, 142 |
| Qualitative peak processing template 112 | sensitivity 196 |
| Quan Flags 338 | Separate ion overlay display 180 |
| Quan mass 339 | Set Chromatogram Reference Sample 109 |
| Quan peak m/z 338 | Set chromatogram reference sample 112 |
| Quan peak response 338 | Set Reference sample 112 |
| Quan peak RT 338 | Shade row when sample is outside of evaluation criteria |
| R^2 threshold 161 | 180 |
| Ranges, compound detection 134 | Show chromatogram on Quantitation Report 180 |
| Raw filename 96 | Showing 337 |
| Recent Files (master methods) 87 | Smoothing |
| Reference compound 120 | Avalon 79 |
| Remove user 35 | Genesis 73, 137 |
| Report concentration 180 | ICIS 76, 140, 143 |
| Report Name, standard reports 176 | Specify default ion ratio ranges 196 |
| Report noise as 75, 138 | Standard type 151 |
| | |

| Status 338 | Polarity parameter |
|---|---|
| Stepoff value 112 | Acquisition List 117 |
| Tailing factor | Compound Datastore 49 |
| Genesis 74, 138 | P-P Threshold, event type 81 |
| ICIS 77, 140 | Precursor Mass parameter |
| Theo Conc | Acquisition List 116–117 |
| Lab Control page 168 | Compound Datastore 49–50 |
| Matrix Spike page 170 | Print parameter, standard reports 176 |
| Method Val page 169 | procedures |
| Surrogate page 167 | |
| Theoretical amount 338 | acquisition mode |
| Time 80 | access real-time display 246 |
| Total response 338 | add samples to the samples list 224 |
| Total rows 337 | assign a specific channel to a sample 229 |
| Trace, Detection 133 | create a batch template 223 |
| Tune file lifetime 181 | export reports to a folder 234 |
| ULOL (Linearity limit) 160 | import samples into the sample list 227 |
| Units 151 | insert samples into the samples list 225 |
| Upper Limit 165 | pause all batches in a queue 253 |
| Use alternate calibration report format 180 | preview a standard report 233 |
| Use an existing raw data file 96 | re-inject a sample from 228 |
| Use as RT Reference 120 | remove a batch from a queue 255 |
| Use autosampler 96 | remove a pending batch from a queue 255 |
| <u> </u> | remove a pending sample from a queued batch 256 |
| Use data dependent scans 198 | remove a single batch from a queue 254 |
| Use Genesis algorithm for qual processing 197 | remove all batches in a queue 254 |
| Use Method Forge 90, 93 | remove all pending batches in a queue 254 |
| Use the default template 96 | remove pending samples from a queued batch 256 |
| Use these libraries 196 | remove samples from the samples list 228 |
| Username 9, 34 | save a to-be-run batch 240 |
| Valley rise (%) 75, 138 | select a previously acquired batch 222 |
| Valley S/N 75, 138 | select a ready-to-acquire batch 221 |
| Value 80 | select a template for a new batch 221 |
| Vial position 96 | specify a calibration batch 238 |
| View width, Detection 72, 131 | specify device states 239 |
| Weighting | specify startup or shutdown methods 237 |
| Calibration 151 | start a new batch 220 |
| Method Template Editor 197 | start an acquisition 240 |
| Window | view output files 243 |
| Acquisition List 117 | configuration mode |
| Compound Datastore 49 | access the Configuration mode 27 |
| Detection 72, 131 | add a compound to the datastore 52 |
| Password parameter | add a confirming ion to a quan peak 54 |
| logon screen 9 | add a quan peak to a compound 53 |
| user administration 34 | add a user 30 |
| password, administrator 8 | change the default drive 40 |
| Path parameter 96 | choose a drive 40 |
| Peak height (%) parameter | copy the folder hierarchy from another drive 43 |
| Genesis 74, 137 | create a new compound datastore 46 |
| ICIS 77, 140 | create projects or subprojects 42 |
| Peak noise factor parameter 77, 140 | delete projects or subprojects 42 |
| Peak S/N cutoff parameter 75, 138 | edit user information 31 |
| Percentage parameter 164 | enable multiplexing 70 |
| Phone number parameter 34 | enable user security 66 |
| Thore number parameter JT | • |

| hide a drive from display 41 | open the Data Review mode 284 |
|---|--|
| import compounds 51 | open the Local Method View 351 |
| import new report types 59 | open the Report View 325 |
| open a compound datastore 46 | print a report 333 |
| open the Application Configuration view 56 | reinject a sample 268 |
| open the Compound Datastore editor 45 | reinject a sample from a previously acquired batch |
| open the Defaults page 68 | 272 |
| open the Project Administration view 39 | remove a manually created peak |
| open the User Administration view 30 | confirming ions 309 |
| open the User Security page 66 | quantification peak 300 |
| refresh the drives display 41 | remove a peak from the peaks pane 315 |
| remove a compound 52 | remove a peak from the qual peak pane 320 |
| remove a confirming ion 54 | remove samples from the samples list 267 |
| remove a user 33 | scroll the samples list 286 |
| remove all empty folders 43 | search for text 335 |
| save a datastore 46 | select a compound 328 |
| save a datastore to a new name 47 | select a report |
| specify common detection parameters 71 | generating reports 330 |
| specify default laboratory and instrument names 69 | viewing reports 326 |
| specify default mass precision and intensity scale 69 | select a sample |
| data review mode | generating reports 331 |
| access the Data Review mode 261 | |
| | viewing reports 328 |
| add samples to the samples list 265 | sort the compounds list 296 |
| change the displayed information for a detected | submit all samples in the batch 278 |
| peaks 309 | submit selected samples 279 |
| change the displayed information for detected peaks | switch between method and manual integration |
| 301, 320 | qual mode 320 |
| change the library entry for a selected peak 323 | quan mode 301 |
| copy a sample in the samples list 267 | temporarily customize the column display |
| create a new batch 264 | Batch view 270 |
| customize the column display 269, 287 | Data Review view 286 |
| display peaks for a specific compound 297, 314 | zoom in on a peak |
| display specific problems with a compound 297 | chromatogram navigation 317 |
| display the internal standard for a compound 298 | confirming ions 309 |
| display the peak information window 310 | ion overlay 312 |
| edit sample values 268 | qual peak 319 |
| edit samples in a batch 272 | quantification peak 299 |
| edit the batch-level output formats 275 | spectra pane 322 |
| edit the sample-level output formats 274 | zoom in on an area |
| enlarge the report text 335 | calibration curve 305 |
| exclude a calibration point 291 | QED spectra 307 |
| export a report 334 | spectra 306 |
| insert samples in the samples list 266 | getting started |
| make a compound active or inactive 290 | choose a node 22 |
| manually add a peak | display a log of instrument errors 23 |
| chromatogram pane 317 | install the NIST library 10 |
| confirming ions 309 | install the QED library 10 |
| qual peak pane 319 | install the TraceFinder application 8 |
| quantification peak 300 | log on to the TraceFinder application 8 |
| manually exclude a calibration point 305 | monitor the instrument status 23 |
| manually integrate a quan or qual ion 300 | start the TraceFinder application 8 |
| modify the peak detection settings 301 | watch the real-time display from the dashboard 24 |
| open a recent batch 272 | 1 / |
| open a saved batch 271 | |
| | |

| method development mode 152 | replace a quan mass 125 |
|--|--|
| access the Method Development mode 85 | resize or reorganize the columns 211 |
| acquire selected samples 212 | save the method template 194 |
| acquire the batch 212 | save the new method 128 |
| add a compound to the list 115 | select compounds from the datastore 102 |
| add a mass as a new compound 126 | set automated background subtraction options 108 |
| add a mass as a new confirming ion 128 | specify a chromatogram reference sample 109 |
| add a mass to the existing quan mass ranges 125 | specify a location for development batch data 208 |
| add a quan peak 126 | specify general information for a master method 106 |
| add a quan peak to an existing compound 146 | specify ion ratio criteria 149 |
| add compounds to the method 100, 122 | specify mass tolerance 110 |
| add confirming ions to an existing compound 147 | specify no maximum concentration 163 |
| add ions to get an accumulated signal 146 | specify peak criteria 189 |
| add sample to the development batch 208 | specify QC levels and concentrations 154 |
| | |
| apply current peak detection settings | specify qualitative peak processing 193 |
| to all peaks in method 135 | specify quantitation flag options 178 |
| to all peaks that use the same detection mode 135 | specify quantitation limits 177 |
| associate a raw data file with the method 99 | specify ranges of ions for detection and integration |
| automatically select compounds to create a new | 132 |
| method 90 | specify standard report types and output formats 173 |
| calculate and report semi-quantitative results 183 | specify the default parameters 182 |
| calibrate the compounds 191 | specify the exact mass window 184 |
| change the compound reference spectrum 125 | specify the Exactive parameters 184 |
| change the quantitation mass used for a quan peak | specify the ion ratio calculation method 184 |
| 144 | specify the maximum concentration as a percentage |
| copy a sample 209 | 163 |
| correct surrogates 179 | specify the maximum concentration as an absolute |
| create a group 171 | value 163 |
| create a method template 188 | specify user interface options 178 |
| create a new instrument method 201 | track the use of the tune file 179 |
| create a new multiplexing instrument method 203 | update confirming ion ratios 144 |
| delete a compound from the list 115 | zoom in the chromatogram or spectrum displays 147 |
| edit the instrument method parameters 107 | Product Mass parameter |
| enter a note for the method 194 | Acquisition List 116–117 |
| enter column values 210 | Compound Datastore 49–50 |
| export SRM data to an XML file 199 | Project Administration view 39 |
| filter the displayed compounds 119, 122 | projects and subprojects, creating 39 |
| identify the peaks 190 | projects and subprojects, creating 39 |
| import a master method 198 | • |
| import an instrument method 206 | Q |
| import an Xcalibur method 97 | QAQC Flags parameter 338 |
| insert samples into the development batch 209 | QAQC role 38 |
| manually select compounds to create a new method | |
| 93 | QC Levels page, Master Method View 154 |
| open a saved master method 104 | QC1-QCn parameter 155 |
| open a saved master method from the Recent Files | QED library, installing 10 |
| list 105 | QED Spectra page 307 |
| | QIon parameter 339 |
| open an instrument method 205 | Qual Mode 284 |
| open the Development Batch view 207 | Qualitative peak processing template parameter 112 |
| open the Instrument View 200 | Quan Flags parameter 338 |
| open the Qual Browser 213 | Quan mass parameter 339 |
| remove all samples from the samples list 211 | |
| remove selected samples from the samples list 211 | Quan Mode 284 |
| replace a confirming ion 128 | Quan peak m/z parameter 338 |

| Quan peak response parameter 338 | Matrix Blank page 163 | |
|---|---|--|
| Quan peak RT parameter 338 | Matrix Spike page 170 | |
| quan report settings, specifying 177, 182 | Method Val page 169 | |
| | QC levels page 155 | |
| R | Solvent Blank page 165 | |
| | Surrogate page 167 | |
| R^2 threshold parameter 161 | | |
| Ranges parameter, compound detection 134 | S | |
| Ratios page, Master Method View 149 | S/N threshold parameter 74, 137 | |
| Raw filename parameter 96 | Sample comment parameter 96 | |
| Real Time Status command 25 | Sample Concentration parameter 339 | |
| Real Time Viewer page, Master Method View 156 | Sample definition view 224 | |
| real-time display 246 | | |
| Recent Files parameter (master methods) 87 | Sample file parameter 337 | |
| Reference compound parameter 120 | sample types, defined 259 | |
| reference spectra 322 | Sample Units parameter 339 | |
| Remove user parameter 35 | Save As Compound Datastore command 47 | |
| Report concentration parameter 180 | Save changes parameter 35 | |
| report formats, specifying 173 | Save Compound Datastore command 46 | |
| Report Name parameter, standard reports 176 | Save Compound Datastore dialog box 47 | |
| Report noise as parameter 75, 138 | Select a custom template parameter 96 | |
| Report selection view 233 | Select a report parameter 337 | |
| Report Title parameter, standard reports 176 | Select batch location parameter 87 | |
| Report View 325 | Select Compounds from CDS parameter 103 | |
| reports | Select Compounds to Add dialog box 103 | |
| Acquisition mode 233 | Selected Reaction Monitoring 55 | |
| configuring 58 | Sensitivity parameter | |
| Data Review mode 325 | Avalon 79 | |
| flags defined 356 | Genesis 73, 137 | |
| listed 5 | ICIS 76, 140, 142 | |
| sample PDFs 357 | sensitivity parameter 196 | |
| viewing landscape in PDF 357 | Separate ion overlay display parameter 180 | |
| Reports page, Master Method View 173 | Set Chromatogram Reference Sample parameter 109 | |
| Reset scaling command 148 | Set chromatogram reference sample parameter 112 | |
| Response parameter 339 | Set Reference sample parameter 112 | |
| Response via parameter | Set this mass as a new quan peak command 148 | |
| Calibration 151 | Set this mass as quan mass command 148 | |
| Method Template Editor 197 | Shade row when sample is outside of evaluation criteria | |
| Retention Time parameter 339 | parameter 180 | |
| RMS parameter 78, 141 | Shoulders On, event type 82 | |
| Role parameter 34 | Show chromatogram on Quantitation Report parameter 180 | |
| RT parameter | Show Internal Standard command 298 | |
| Acquisition List 117 | Showing parameter 337 | |
| Calibration 151 | SIM experiment type 55 | |
| Calibration levels 153 | Single Ion Monitoring 55 | |
| Calibration page 161 | Smoothing parameter | |
| Chk Std page 162 | Avalon 79 | |
| Compound Datastore 49 | Genesis 73, 137 | |
| Identification 119 | ICIS 76, 140, 143 | |
| ISTD page 166 | Solvent Blank page, Master Method View 165 | |
| Lab Control page 168 | Specify default ion ratio ranges parameter 196 | |
| Limits page 160 | | |

| Spectra page 306 | Use an existing raw data file parameter 96 |
|---|---|
| Spectrum page, Master Method View 144 | Use as RT Reference parameter 120 |
| SRM data, exporting 199 | Use autosampler parameter 96 |
| SRM experiment type 55 | Use data dependent scans parameter 198 |
| standard reports | Use Genesis algorithm for qual processing parameter 197 |
| listed 6 | Use Method Forge parameter 90, 93 |
| standard reports, list of 5 | Use the default template parameter 96 |
| Standard type parameter 151 | Use these libraries parameter 196 |
| Start Threshold, event type 81–82 | user accounts, creating 30 |
| starting TraceFinder 8 | User Administration view 30 |
| Status parameter 338 | user roles and permissions 36 |
| Stepoff value parameter 112 | user roles, defined 36 |
| Supervisor role 37 | user security, enabling 66 |
| supported file types, defined 2 | Username parameter 9, 34 |
| supported software and hardware viii | commine parameter 7, 31 |
| Surrogate page, Master Method View 167 | V |
| system requirements viii | |
| 1 | Valley rise (%) parameter 75, 138 |
| Т | Valley S/N parameter 75, 138 |
| | Value parameter 80 |
| Tailing factor parameter | Vial position parameter 96 |
| Genesis 74, 138 | View width parameter, Detection 72, 131 |
| ICIS 77, 140 | views |
| Tangent Skim, event type 82 | Batch 262 |
| target screening reports, list of 6 | Batch/Template Selection 220 |
| Technician role 37 | Compound Datastore 48 |
| templates | Data Review 284 |
| batch 223 | Development Batch 207 |
| method 188 | Finish 237 |
| Tension, event type 82 | Instrument 200 |
| Theo Conc parameter | Local Method 351 Method 89 |
| Lab Control page 168 | Project Administration 39 |
| Matrix Spike page 170 | Report selection 233 |
| Method Val page 169 | Report View 325 |
| Surrogate page 167 | Sample definition 224 |
| Theoretical amount parameter 338 | User Administration 30 |
| Thermo Xcalibur Instrument Setup dialog box 107 | |
| Time parameter 80 | W |
| Total response parameter 338 | |
| Total rows parameter 337 | Weighting parameter |
| Trace parameter, Detection 133 | Calibration 151 |
| Tune file lifetime parameter 181 | Method Template Editor 197 |
| | Window parameter |
| U | Acquisition List 117 |
| ULOL (Linearity limit) parameter 160 | Compound Datastore 49 |
| Units parameter 151 | Detection 72, 131 |
| Unknown sample type, defined 259 | workflow |
| Unknown/TIC sample type, defined 259 | acquisition mode 218 |
| Update confirming ion ratios with this spectrum command | general 4 |
| 148 | V |
| Upper Limit parameter 165 | X |
| Use alternate calibration report format parameter 180 | XIC experiment type 55 |

Index: X