

Thermo Xcalibur

Quan Browser

User Guide

Software Version 3.1

XCALI-97615 Revision A August 2014



Thermo
SCIENTIFIC

© 2014 Thermo Fisher Scientific Inc. All rights reserved.

Xcalibur is a registered trademark of Thermo Fisher Scientific Inc. in the United States.

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

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, August 2014

Software version: Xcalibur 3.1 and later

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

Contents

| | | |
|------------------|---|------------|
| | Preface | vii |
| | Related Documentation | vii |
| | Safety and Special Notices | viii |
| | Contacting Us | ix |
| Chapter 1 | Overview of Quantitative Analysis | 1 |
| | Acquiring and Quantitatively Processing Data | 2 |
| | Integrating and Identifying Chromatographic Peaks | 4 |
| | Using External Standards | 6 |
| | Using Internal Standards | 8 |
| Chapter 2 | Quan Browser Overview | 11 |
| | Understanding How Quan Browser Works | 11 |
| | Calibration Replicates | 11 |
| | Named Calibration File | 12 |
| | Brackets and Groups for Sequences | 12 |
| | Unbracketed Sequence | 12 |
| | Open Bracket Sequence | 14 |
| | Nonoverlapping Bracket Sequence | 14 |
| | Overlapping Bracket Sequence | 14 |
| | Starting Quan Browser | 15 |
| | Quan Browser Window Overview | 17 |
| | Component List | 17 |
| | Results Grid | 18 |
| | Chromatogram Plot and Companion Views | 19 |
| | Saving Changes Made in Quan Browser | 21 |
| Chapter 3 | Working with the Quan Browser Results Grid | 25 |
| | Reviewing and Reworking the Results Grid | 25 |
| | Reviewing the Peak Status Column | 27 |
| | Making Changes in the Results Grid | 27 |
| | Adding and Removing Samples | 28 |
| | Hiding or Displaying Columns | 31 |
| | Changing the Sort Order | 32 |

| | | |
|-------------------|--|-----------|
| Chapter 4 | Working with Peak Identification and Detection | 33 |
| | Viewing and Changing Peak Identification, Detection, and Integration | 33 |
| | Setting the Display Options for the Chromatogram and Spectrum Plot Views | 35 |
| | Viewing Peak Information | 35 |
| | Using the User Identification Settings Dialog Box | 37 |
| | Setting the Identification Values | 38 |
| | Setting the Detection Parameters | 40 |
| | Setting the Integration Parameters | 41 |
| | Setting the Advanced Parameters for the Genesis and ICIS Algorithms | 43 |
| | Setting the Flags Values | 45 |
| | Integrating Chromatogram Peaks Manually | 46 |
| | Using the Spectrum Plot Companion View | 48 |
| | Reviewing Spectrum Search Results for GC/MS Data | 50 |
| Chapter 5 | Working with the Calibration Settings | 53 |
| | Reviewing and Reworking the Calibration Settings in Quan Browser | 53 |
| | Reviewing and Reworking the Calibration Curve Settings | 55 |
| | Changing the Isotope Percentage Values | 57 |
| | Excluding Data Points from the Calibration Curve | 59 |
| Chapter 6 | Viewing Results and Generating Reports | 65 |
| | Reviewing the Results of the Quantitative Analysis | 65 |
| | Generating Reports | 68 |
| Appendix A | Quan Browser Window | 73 |
| | Quan Browser Window | 74 |
| | Quan Browser Title Bar | 75 |
| | Quan Browser Menu Bar | 76 |
| | File Menu – Quan Browser | 76 |
| | View Menu – Quan Browser | 78 |
| | Zoom Menu – Quan Browser | 79 |
| | Options Menu – Quan Browser | 79 |
| | GoTo Menu – Quan Browser | 80 |
| | Help Menu – Quan Browser | 80 |
| | Quan Browser Toolbar | 81 |
| | Quan Browser Results Grid | 84 |
| | Quan Browser Component List | 88 |
| | Quan Browser Chromatogram Plot View | 88 |
| | Quan Browser Spectrum Plot View | 90 |
| | Quan Browser Calibration Curve View | 92 |

| | |
|---|------------|
| Quan Browser Dialog Boxes | 94 |
| Add Sample Dialog Box | 95 |
| Bracket/Group In Use List | 96 |
| Cal Exclusion List Dialog Box | 97 |
| Calibration Settings Dialog Box | 98 |
| Curve Page – Calibration Settings Dialog Box | 98 |
| Flags Page – Calibration Settings Dialog Box | 100 |
| Isotope% Page – Calibration Settings Dialog Box | 101 |
| Levels Page – Calibration Settings Dialog Box | 104 |
| Type Page – Calibration Settings Dialog Box | 105 |
| Display Options Dialog Box in Quan Browser | 107 |
| Masses Dialog Box | 107 |
| Peak Information Dialog Box | 108 |
| No Peak Page – Peak Information Dialog Box | 109 |
| Info Page – Peak Information Dialog Box | 110 |
| Flags Page – Peak Information Dialog Box | 111 |
| More Flags Page – Peak Information Dialog Box | 114 |
| Chro Page – Peak Information Dialog Box | 115 |
| Info or More Info Page – Peak Information Dialog Box | 116 |
| More Info Page – Peak Information Dialog Box | 117 |
| Spectrum Page – Peak Information Dialog Box | 118 |
| Suitability Page – Peak Information Dialog Box | 118 |
| Quantitation Results Sorting Order Dialog Box | 119 |
| Reports Dialog Box | 121 |
| Result List Column Hiding Dialog Box | 126 |
| Select Level Dialog Box | 127 |
| Select Report Samples Dialog Box | 128 |
| User Identification Settings Dialog Box | 129 |
| Identification Page – User Identification Settings Dialog Box | 130 |
| Detection Page – User Identification Settings | 134 |
| Integration Page – User Identification Settings Dialog Box | 139 |
| Genesis Integration Page Parameters | 139 |
| ICIS Integration Page Parameters | 141 |
| Avalon Integration Page Parameters | 142 |
| Advanced Page – User Identification Settings Dialog Box | 145 |
| Genesis Advanced Page Parameters | 145 |
| ICIS Advanced Page Parameters | 147 |
| Flags Page – User Identification Settings Dialog Box | 148 |
| View Sample Types Dialog Box | 149 |
| Index | 151 |

Preface

This guide describes how to use the Thermo Xcalibur™ Quan Browser window to review and rework the results of a quantitative analysis.

Contents

- [Related Documentation](#)
- [Safety and Special Notices](#)
- [Contacting Us](#)

To provide us with comments about this document, please click the link below. Thank you in advance for your help.



Before using this guide, read your instrument's *Getting Started Guide* and the *Xcalibur Data Acquisition and Processing User Guide* to become familiar with the basic features of the Xcalibur data system such as the Home Page, Instrument Setup, and Processing Setup windows.

Related Documentation

Thermo Fisher Scientific provides the following documentation for the Xcalibur data system:

- *Xcalibur Quantitative Analysis Getting Started Guide*
- *Xcalibur Data Acquisition and Processing User Guide*
- *Xcalibur Quan Browser User Guide*
- *Xcalibur Qual Browser User Guide*
- *Xcalibur Library Browser User Guide*
- *XReport User Guide*
- Help from within the data system

❖ **To access the manuals from the data system computer**

- From the Windows™ Start menu, choose **All Programs (or Programs) > Thermo Xcalibur > Manuals**.
- From the Home Page – Roadmap view of the Thermo Xcalibur data system, choose **Help > Manuals** from the menu bar.

Thermo Fisher Scientific provides the manuals as PDF files.

Safety and Special Notices

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

Safety and special notices include the following:

IMPORTANT Highlights information necessary to avoid damage to software, loss of data, invalid test results, or information critical for optimal performance of the system.

Note Highlights information of general interest.

Tip Highlights helpful information that can make a task easier.

Contacting Us

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

| For Thermo Scientific™ products | Access by phone, fax, email, or website |
|---|--|
| Technical Support | <p>(U.S.) Phone: 1 (800) 532-4752 Fax: 1 (561) 688-8736</p> <p>Email: us.techsupport.analyze@thermofisher.com</p> <p>Web—for product support, technical documentation, and knowledge bases: www.thermoscientific.com/support</p> |
| Customer Service (Sales and service) | <p>(U.S.) Phone: 1 (800) 532-4752 Fax: 1 (561) 688-8731</p> <p>Email: us.customer-support.analyze@thermofisher.com</p> <p>Web—for product information: www.thermoscientific.com/lc-ms</p> <p>Web—for customizing your service request:</p> <ol style="list-style-type: none"> 1. From any Products & Services web page, click Contact Us. 2. In the Contact Us box, complete the information requested, scroll to the bottom, and click Send. |
| User Documentation | <p>Web—for downloading documents: mssupport.thermo.com</p> <ol style="list-style-type: none"> 1. On the Terms and Conditions web page, click I Agree. 2. In the left pane, click Customer Manuals. 3. To locate the document, click Search and enter your search criteria. For Document Type, select Manual. <p>Email—to send feedback directly to Technical Publications: techpubs-lcms@thermofisher.com</p> <p>Web—to complete a survey about this Thermo Scientific document: www.surveymonkey.com/s/PQM6P62</p> |

Overview of Quantitative Analysis

The Xcalibur data system is a complete quantitative and qualitative analysis software package that you can use to acquire data specifically for analytes of interest, to perform confirmatory library searches, and to determine the concentration of analytes in samples. The Xcalibur data system interfaces with the XReport reporting package to print individual sample reports and sequence summary reports for analyses. For more information on XReport, refer to the *XReport User Guide*.

Quan Browser is a powerful and versatile utility for reviewing and reworking the following:

- Component peak identification and integration criteria
- Standards, QCs, blanks, and unknowns
- Calibration curves for quantitation standards

After making any changes, save the new results with an audit trail describing the reason for the change.

Quan Browser incorporates a calibration curve display, peak integration, and results view where you can do the following:

- Process quantitation sequences
- Interactively edit processing parameters and audit the changes
- Create new files that keep track of processing results for individual raw files and include a copy of the method used to generate the results

Result files changed using Quan Browser do not affect the original processing method.

This chapter describes some of the basic principles and terminology of quantitation, and provides a brief overview of quantitation with the Xcalibur data system.

Contents

- [Acquiring and Quantitatively Processing Data](#)
- [Integrating and Identifying Chromatographic Peaks](#)

Acquiring and Quantitatively Processing Data

Quantitative analysis is the process of measuring the amount of a particular component in a sample. With the Xcalibur data system, quantitative analysis usually involves the following steps:

Note The order of some of these steps is not rigid. For example, you can acquire and process a set of data files using a sequence that contains both an instrument method and a processing method, and you can print reports without previewing them first. For more information, refer to the *Xcalibur Data Acquisition and Processing User Guides*.

1. Create an instrument method.

The Xcalibur data system uses an instrument method to store a specific set of parameters used to operate the autosampler, LC pump or MS pump, gas chromatograph, mass spectrometer, PDA detector, and so on.

For more information about creating an instrument method, refer to your hardware documentation.

2. Create an acquisition sequence.

An acquisition sequence identifies the position of the samples in an autosampler tray (if appropriate), the instrument method used to control the HPLC, GC/MS, or LC/MS instrument, and the directory and file names for the acquired data files.

For more information about creating an acquisition sequence, refer to the *Xcalibur Data Acquisition and Processing User Guide*.

3. Run the sequence to acquire the raw data files.

Run either one sample or a series of samples from the current sequence.

For more information about running samples, refer to the *Xcalibur Data Acquisition and Processing User Guide*.

4. Create a processing method.

Create processing methods in the Processing Setup view. The data system uses a processing method to identify, detect, and integrate components in a chromatogram, generate calibration curves, quantify unknowns, and produce reports. The application contains several built-in report templates. Report templates have an (.xrt) file extension.

For more information about creating a processing method, refer to the *Xcalibur Data Acquisition and Processing User Guide*.

5. Create a processing sequence by adding the processing method to the original acquisition sequence.

A processing sequence contains a processing method, consists of a list of sample data files, and includes information on sample type and calibration or QC level.

6. Process a representative raw file or the entire sequence with the processing method by using the Batch Reprocess feature in the sequence Setup view.

Processing a raw file produces a result file. Result files have an (.rst) file extension. For instructions on batch processing a sequence, refer to the *Xcalibur Data Acquisition and Processing User Guide*.

Tip The report templates provided with the Xcalibur software are generic and might not produce the results you expect. Preview a report in XReport before printing reports for an entire sequence. For information about creating and changing reports, refer to the *XReport User Guide*.

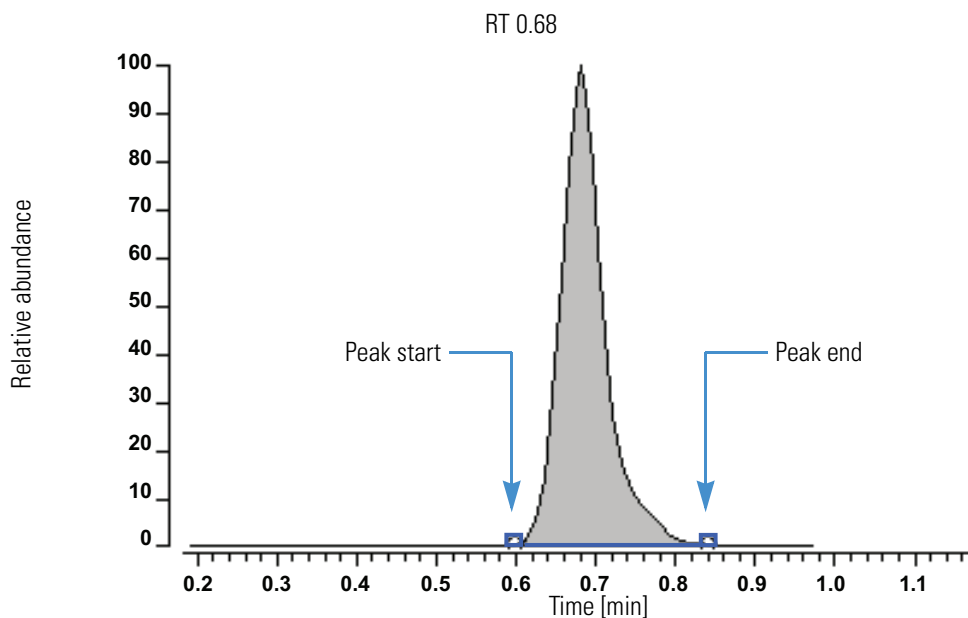
7. Review the processed data files. After the Xcalibur data system processes the raw data files, open the processed sequence in Quan Browser and evaluate the peak detection settings, the integration settings, and the calibration curve for each named component. As you evaluate the results of the processing method, you can modify some of its parameters in Quan Browser. If the processing method contains a report template, you can print reports from Quan Browser.
8. Preview a report for a representative data file from the XReport reporting package.

To produce customized reports, open a representative result file (RST) in the XReport reporting package and create a report template.
9. Once you are satisfied with the way a report displays your data, add the report to the processing method, if you have not already done so, and batch process the sequence to generate printed reports.

Integrating and Identifying Chromatographic Peaks

The Xcalibur data system integrates chromatograms to separate the chromatographic peaks from the baseline noise, identify the beginning and end of each peak, identify the peak maxima, and calculate the area or height of each peak (see [Figure 1](#)). The data system uses one of the following three algorithms to detect and integrate the peaks in chromatograms: Genesis, ICIS, or Avalon. To integrate mass chromatograms, use either the Genesis (designed for legacy Xcalibur 1.0 studies) or the ICIS integration algorithm. To integrate UV/Vis and analog chromatograms, use the Avalon integration algorithm.

Figure 1. Integrated chromatographic peak, showing peak start and peak end markers

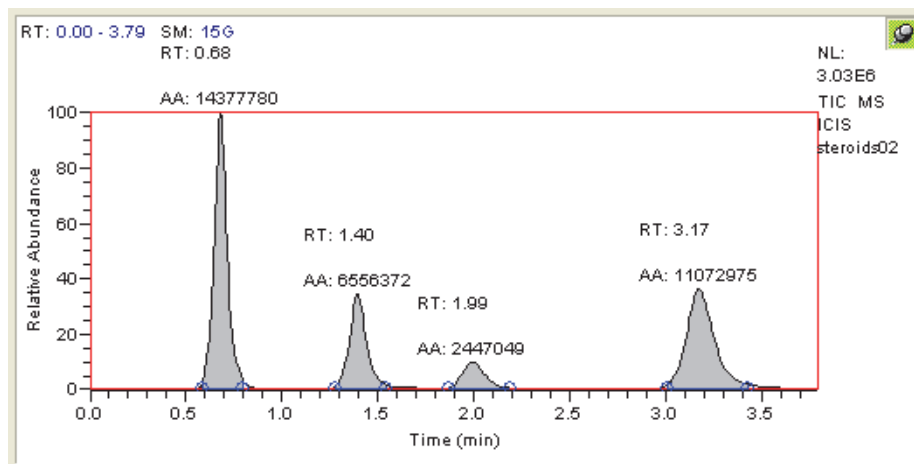


For LC data, the Xcalibur data system identifies peaks based on their retention times. The retention time of a peak is the time that elapses between the injection of the sample and the detection of the peak maxima. For GC data, the application can identify peaks based on either their retention times or their mass spectra.

During a sequence run, the retention times of chromatographic peaks can vary slightly. As a result, you must enter an appropriate retention time window for each peak, in addition to its expected retention time. A retention time window is a time range that brackets the discrete retention time setting. The appropriate retention time window for a chromatographic peak depends on several factors, including the width of the chromatographic peak and the specificity of the chromatographic method. Due to band broadening as the sample travels through the column, highly retained compounds produce wider chromatographic peaks. So in general, use a wider retention time window for late eluting compounds than for early eluting compounds. [Figure 2](#) shows the effect of retention time on peak width. The

chromatogram shown in Figure 2 contains four peaks. The retention times and widths of these peaks are listed in the following table. In this example, hydrocortisone, which elutes at 0.68 min, has a peak width of 0.2 min; whereas, progesterone, which elutes at 3.17 min, has a peak width of 0.6 min.

Figure 2. Integrated total ion current chromatogram (TIC) for steroids02.raw



| Compound | Retention time (min) | Baseline peak width (min) |
|---------------------|----------------------|---------------------------|
| hydrocortisone | 0.68 | 0.2 |
| deoxycorticosterone | 1.40 | 0.3 |
| methyltestosterone | 1.99 | 0.4 |
| progesterone | 3.17 | 0.6 |

Because the Xcalibur data system might detect more than one chromatographic peak within the specified retention time window, identify the target compound as either the highest peak in a chromatogram or the closest peak to the expected retention time. Use the Genesis and ICIS integration algorithms (used for mass spectral data) to rule out peaks below a specified signal-to-noise ratio.

Quantitative analysis of samples containing unknown amounts of the target component is achieved by first calculating the peak area or height and then computing and applying the appropriate response to the equation derived from the calibration curve. This process provides an estimate of the amount of the unknown component. The precision of the measurement depends on the quality and, to a lesser extent, the quantity of the calibration data.

The detection limit of the quantitation method is the lowest concentration of analyte in a sample that can be detected but not necessarily calculated as an exact value. The lower and upper quantitation limits are the lowest and highest concentrations of analytes in a sample that can be measured with an acceptable level of accuracy and precision, respectively. In an analytical method, the highest concentration calibration standard defines the upper

quantitation limit. The quantitation range is the range of concentration between the lower and upper quantitation limits (including these limits) that can be reliably quantified time after time with acceptable levels of accuracy and precision through the use of a concentration-response relationship.

There are two basic quantitation techniques:

- External standard (ESTD) quantitation
- Internal standard (ISTD) quantitation

The chosen method determines the calculation method, both for the generation of the calibration curves and for subsequent quantitation.

For information about using the two basic quantitation techniques, see these topics:

- [Using External Standards](#)
- [Using Internal Standards](#)

Using External Standards

An external standard (ESTD) is a separate sample that contains a known amount of the target compound. To perform an ESTD calibration, prepare a set of standard solutions containing a known amount of the target compounds. After you inject these solutions, the data system analyzes the resulting chromatograms and constructs a calibration curve for each target compound by plotting the magnitude of the detector's response as a function of the amount of the target compound according to the following equation:

$$\text{Response}_{\text{cal}} = f(\text{Amount}_{\text{cal}})$$

Where:

f = curve type

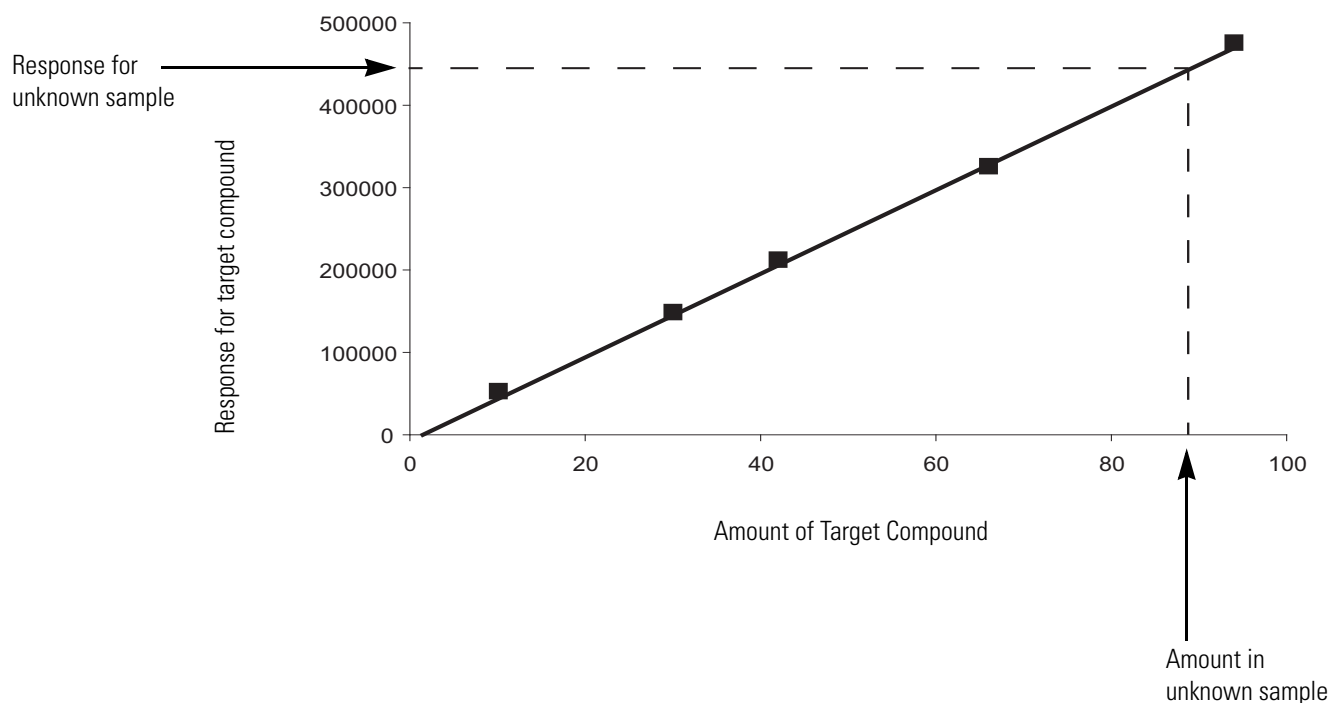
$\text{Amount}_{\text{cal}}$ = amount of calibration standard

$\text{Response}_{\text{cal}}$ = response for calibration standard

The data system determines the amount of the target compounds in each unknown by comparing the magnitudes of their responses to the calibration curves (see [Figure 3](#)).

Use ESTDs if all compounds of interest can be assayed by using a single set of external standards. This approach offers time- and cost-effective quantitation for applications using high precision autosamplers and traditional UV/Vis detectors. However, for some types of analyses, this method cannot achieve the highest level of precision and accuracy. Depending on the instrumentation, variations in analyte and solution stability, injection reproducibility, and matrix interference can lead to lower precision levels in the external standard method than in the internal standard method.

Figure 3. Calibration curve generated by using an external standard



In general, the external standard calibration is an effective quantitation technique; however, if one or more of the following problems exist, consider using the internal standard calibration technique instead.

- Lack of injection reproducibility
- Changes in analyte solution volume
- Matrix and coeluter interference (both suppression and enhancement)
- System instability
- Variations in the source conditions

Using Internal Standards

An internal standard (ISTD) is a component that is added to a sample to act as a response reference for one or more non-ISTD components in the sample. The concentration or amount of an ISTD in any standard or unknown sample typically remains constant.

Because quantitative mass spectrometric analysis usually involves multiple steps, the total error in the analysis results from the accumulation of errors at each step. In general, sample handling errors account for a larger fraction of the total error than detector errors do. Fortunately, the internal standard method can reduce both sources of error. For example, internal standards can correct for variations in a component's peak area that are caused by the following:

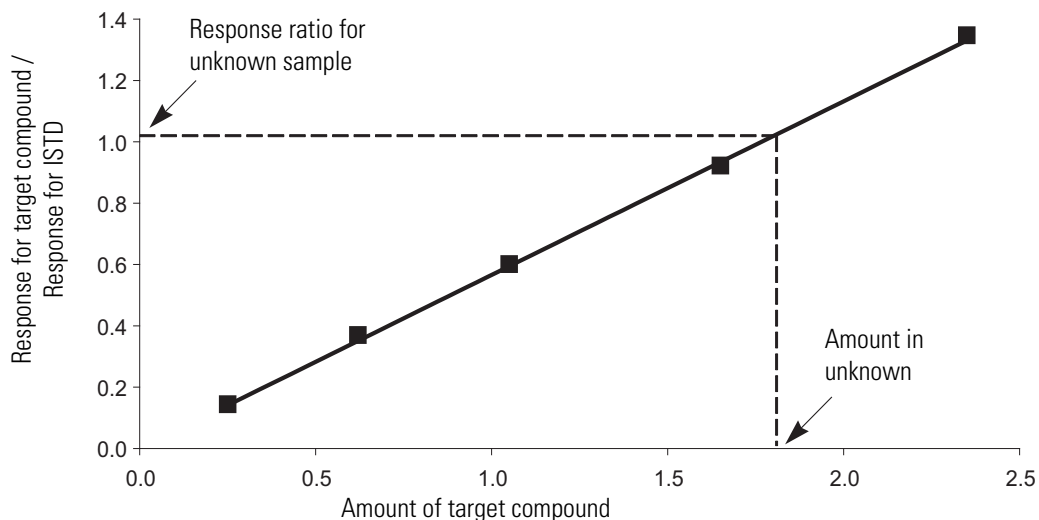
- Lack of injection reproducibility
- Changes in analyte solution volume
- Matrix and coeluter interference (both suppression and enhancement)
- System instability
- Variations in the source conditions

For maximum precision, add the ISTD component as early as possible to the start of the sample workup, particularly in those quantitative methods that require sample manipulations such as extraction, cleanup, and dilution. Since the ISTD and non-ISTD components are analyzed together and since the ISTD is known and quantifiable, the internal standard quantitation approach corrects for injection and other sample handling errors. The ISTD must behave chemically in an identical or similar manner to the target compound through the extraction, cleanup, and analytical processes.

You can also add the ISTD component as the last step of sample preparation prior to the sample's use to compensate for fluctuations in the reproducibility of the sample injection.

For the internal standard calibration method, the Xcalibur data system constructs a calibration curve from a set of standard solutions that contain a range of concentrations for the target compounds and a fixed concentration for each internal standard compound. For each target compound, the data system plots the detector response ratios for the target compound and its associated internal standard compound as a function of the corresponding target compound concentration. The data system then determines the concentration of the target compounds in the unknown samples by using the calibration curve. [Figure 4](#) shows the interpolation of the target compound amount in an unknown from calibration curve constructed using the internal standard method.

Figure 4. Calibration curve generated by using the internal standard method



The calibration curve is determined by the following equation:

$$(\text{Response Ratio}_{\text{TargetCal/ISTD}}) = f(\text{Amount}_{\text{TargetCal}})$$

Where:

$\text{Amount}_{\text{TargetCal}}$ = Amount of target compound in the calibration standards

$\text{Response Ratio}_{\text{TargetCal/ISTD}}$ = ratio of the responses of the target compound to the internal standard compound in the calibration standard

f = equation of the calibration curve according to the selected fit type

Ideally, an ISTD is closely related to the target component in terms of its physical and chemical properties. If the ISTD is used only to compensate for injection reproducibility or changes in the analyte solution volume, it must possess a similar retention (k) to the target component, but it does not need to be chemically similar to the target component. It must be pure, not present in the sample, and inert towards the components of the sample. ISTD components are typically analogs, homologues, or isomers of the target non-ISTD component. An ideal ISTD is a structural or isotopically-labeled analog of one of the target components. Stable isotope-labeled ISTDs act almost identically to the analyte throughout sample manipulation and with regard to ionization tendencies and fragmentation. Internal standards labeled with two or more deuterium (D) atoms are frequently used for LC/MS.

There can be any number of ISTD components in a sample, but each non-ISTD component can be calibrated against only one ISTD component.

Quan Browser Overview

This chapter describes how to use Quan Browser to review processed quantitation sequences. It explains the properties and uses of each component within the Quan Browser window. The *Xcalibur Qual Browser User Guide* describes how to use the Qual Browser window to review raw files. The *Xcalibur Library Browser User Guide* describes how to create user libraries of mass spectral data and run searches of mass spectral libraries.

Contents

- [Understanding How Quan Browser Works](#)
- [Starting Quan Browser](#)
- [Quan Browser Window Overview](#)
- [Saving Changes Made in Quan Browser](#)

Understanding How Quan Browser Works

Using the Quan Browser application, you can step through a sequence of processed samples and review the results for each component in each sample.

The Xcalibur data system has the following quantitation features:

- [Calibration Replicates](#)
- [Named Calibration File](#)
- [Brackets and Groups for Sequences](#)

Calibration Replicates

Calibration replicates are multiple injections of the calibration mixture at the same calibration level or amount. These standard samples all contain the same amount of target compound, so they correspond to the same calibration level. Choose replicates to include or exclude from the calibration curve by using the calibration curve view.

Named Calibration File

After creating a sequence with the Bracket Type set to None, specify a calibration file name in the Calibration File box. Although in theory a different calibration file name for every sample is possible, in practice only one name per sequence is common.

Named calibration files are not available with bracketed sequences.

Brackets and Groups for Sequences

This section describes different bracket types and when to use them to get a specific result.

- [Unbracketed Sequence](#)
- [Open Bracket Sequence](#)
- [Nonoverlapping Bracket Sequence](#)
- [Overlapping Bracket Sequence](#)

To set up sequence bracketing for the None, Non-Overlapped, and Overlapped bracket types, you must use the New Sequence Template dialog box. When you manually set up a sequence using the sequence table in the Sequence Setup view, you can only create an open bracket sequence. You cannot use a calibration file or add more than one processing method to an open bracket sequence. For information about using the New Sequence Template dialog box to set up a sequence, refer to the *Xcalibur Data Acquisition and Processing User Guide*.

Unbracketed Sequence

The Xcalibur data system processes an unbracketed sequence (None bracket type) using a procedure known as the continuing calibration method. Each time the application processes an unbracketed sequence, it creates or updates the calibration files named in the sequence.

Select this process and avoid using Std Clear (Standard Clear) to add replicate data incrementally to a calibration file without discarding the existing replicate data.

The Quan Browser application breaks down unbracketed sequences into logical groups that are somewhat analogous to brackets. It does this by first ordering the samples chronologically with respect to acquisition date and time. It then examines the sequence and starts a new group whenever it encounters a standard. The group ends at the nonstandard sample that immediately precedes the next standard found.

The first group always starts with the first sample, even if it is not a standard. The last group always ends with the last sample. Further, a Std Clear always starts a new group, even if no intervening nonstandard sample has been found following one or more Std Updates.

The Xcalibur data system forms additional logical groups if different named calibration files have been specified in the Cal File entries of the sequence. Each cal file entry causes a new group to be formed. Because using multiple-named calibration files is not typical, their use is not considered any further in this document, but should be deducible from the discussions on groups.

As the Quan Browser application processes each group, it quantifies samples against the current calibration curve. The application processes each standard that it encounters by either replacing (sample type set to Std Clear) or adding to (sample type set to Std Update) the calibration replicate list, generating a new calibration curve.

Quan Browser processing closely emulates that of batch processing (either batch processing directly after acquisition or, subsequently, as a batch process operation). If the Quan Browser application cannot find or open a specified calibration file, the Xcalibur data system displays this message in the Calibration File edit box:

Cal File Unavailable – Using Embedded Calibration

The Quan Browser application takes replicate data from the data stored in the result file. In most cases this data is identical to the data contained within the original calibration file.

Once the Quan Browser application has set up the groups, they are independent and are effectively treated as brackets. In other words, changes in one group do not affect any other group, unlike in batch processing where subsequent groups might well be affected.

The following list illustrates the procedure (for a single-named calibration file):

| | | |
|-----------|------------|-------------------|
| Sample 1 | Unknown | Group 1 start |
| Sample 2 | Unknown | Group 1 end |
| Sample 3 | Std Clear | Group 2 start |
| Sample 4 | Unknown | Group 2 end |
| Sample 5 | Std Update | Group 3 start |
| Sample 6 | Std Update | |
| Sample 7 | Unknown | |
| Sample 8 | Unknown | |
| Sample 9 | Blank | |
| Sample 10 | QC | Group 3 end |
| Sample 11 | Std Update | Group 4 start/end |
| Sample 12 | Std Clear | Group 5 start |
| Sample 13 | Blank | Group 5 end |

Open Bracket Sequence

For an open bracket sequence (Open bracket type), the Quan Browser application creates a replicate list directly from all standard samples in the sequence without using any calibration data embedded in result files.

When you open a single result file in the Quan Browser application, the application treats it as a sequence with only one entry and lists the sample type as Unknown. To show the calibration curve used to quantitate the sample, the application creates the replicate list from the embedded information.

Nonoverlapping Bracket Sequence

For a non-overlapping bracket sequence (Non-Overlapping bracket type), the Quan Browser application creates a separate replicate list for each bracket. The application creates each replicate list directly from all standard samples in the bracket without using any calibration data embedded in result files.

Overlapping Bracket Sequence

For an overlapping bracket sequence (Overlapping bracket type), the Quan Browser application creates a separate replicate list for each bracket. The application creates each replicate list directly from all standard samples in the bracket without using any calibration data embedded in result files.

Exceptions occur for shared standard samples between brackets. When a standard that is shared undergoes a change, that change is reflected in all brackets that contain that standard. When a shared standard sample is deleted, the application deletes the standard sample in all brackets that contain that standard sample and adjusts the replicate lists for all brackets.

When you add a sample as a standard to any bracket, the application adds it to the replicate list automatically. To add a standard sample as a shared standard sample, you must add it separately to each bracket.



The exclusion status of the replicates is independent for each bracket. Even shared samples might be excluded in one bracket but not in another. This is the only exception to a shared sample having identical settings.

Starting Quan Browser

Use the Quan Browser window to review the results of a quantitative analysis.

❖ To start the Quan Browser application

1. Do one of the following:

- On the Home Page – Roadmap view, click the **Quan Browser** icon, .
- From the Instrument Setup window, click the **Home Page** icon, , in the toolbar. Then, click the **Quan Browser** icon.
- From the Home Page or Processing Setup windows, choose **GoTo > Quan Browser**.

At startup, Quan Browser displays the Open dialog box so you can select an existing file. If you do not want to select a file, click **Cancel** in the Open dialog box to close Quan Browser.

The Quan Browser application supports these file types:

- Sequence files (SLD)
- Result files (RST)
- Quan Browser files (XQN)

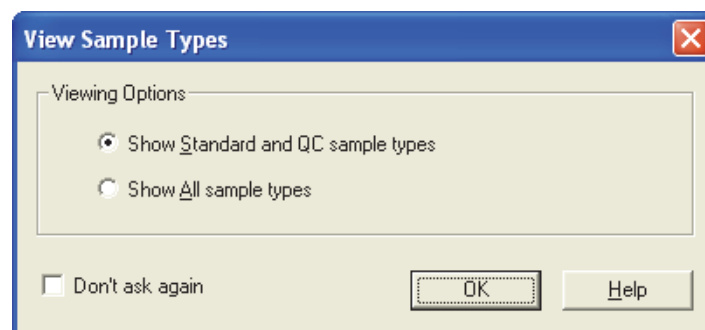
Quan Browser handles result files as single entry sequences.

2. Select a file in the Open dialog box.

When you select a sequence file, the application checks that all the associated raw and result files are available. When it encounters a problem with the sequence file, the application provides information about the likely cause in a warning dialog box and prompts you to exit the application or select a different file.

After the application verifies that the files exist and can be opened, the View Sample Types dialog box opens (see [Figure 5](#)).

Figure 5. View Sample Types dialog box



The two options provided in the View Sample Types dialog box determine how the results grid is configured at startup.

3. Select one of these options:

- To display only Standards and QCs in the Quan Browser results grid, select the **Show Standard and QC sample types** option. Blanks and Unknowns do not appear in the grid. Click either the **Standards** or **QCs** tab.
- To display Standards, QCs, Blanks, and Unknowns in the Quan Browser Grid view, select the **Show All sample types** option. Click one of the following tabs: **All**, **Standards**, **QCs**, **Blanks**, or **Unknowns**.

The View Sample Types dialog box includes a Don't ask again check box. When you select this check box, the dialog box is not displayed when you start subsequent sessions in Quan Browser and the current selection becomes the default.

Note To make this and all other Don't Ask Again-type dialog boxes active, choose **Options > Enable Warnings**.

4. To start the session, click **OK**. Quan Browser loads the specified sequence or file and configures the results grid using your selected viewing option.

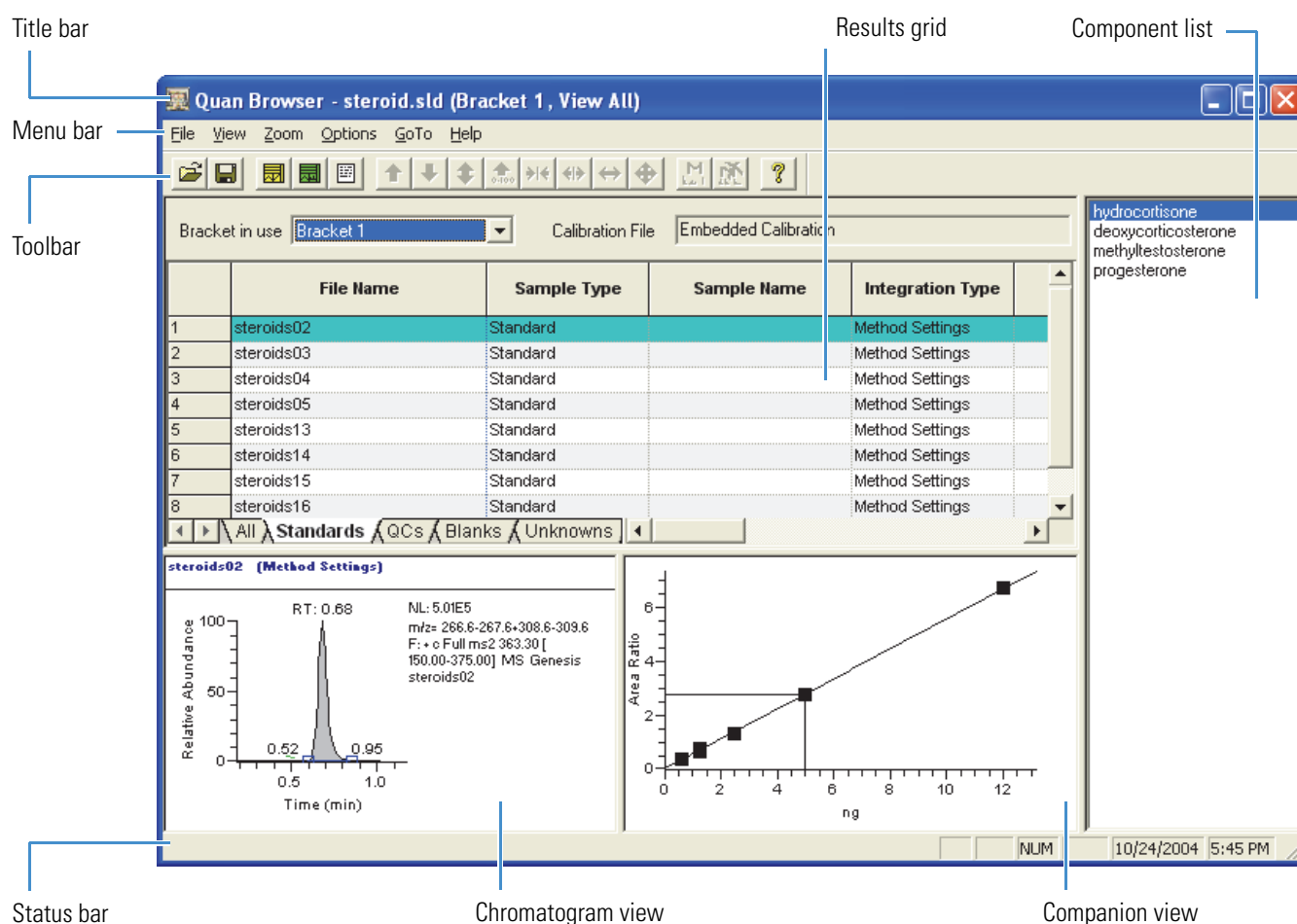
Quan Browser Window Overview

The Quan Browser window has the following features (see [Figure 6](#)):

- [Component List](#)
- [Results Grid](#)
- [Chromatogram Plot and Companion Views](#)

For more information about the Quan Browser window, see [Appendix A, “Quan Browser Window.”](#)

Figure 6. Quan Browser window



Component List

The component list displays all the components within the current bracket sorted by retention time. To update the chromatogram plot view and the companion view with data for a specific component, click the component name. For more information about the component list, see [“Quan Browser Component List”](#) on [page 88](#).

Results Grid

The results grid is made up of the Bracket in Use list, the Calibration File box, and a sequence table. Each row of the sequence table defines a result file and associated parameters. Use the tabs at the bottom of the window to display all of the sequence samples or a subset of the sequence samples.

The results grid is made up of these key areas:

- [Bracket/Group in Use](#)
- [Calibration File](#)
- [Results Grid](#)

Bracket/Group in Use

For bracketed sequences, the Brackets in Use list shows the available brackets in sequential order. The data system selects the first bracket in the list when the file is first loaded into Quan Browser and displays the samples within this bracket in the results grid.

When you load an unbracketed sequence, the samples are broken into logical groups. The Groups in Use list shows the available groups.

Selecting a new bracket or group from the list refills the results grid with the samples from the selected bracket or group. The application updates all the other Views and dialog boxes automatically.

For more information, see [“Brackets and Groups for Sequences”](#) on page 12 and [“Bracket/Group In Use List”](#) on page 96.

Calibration File

This read-only box shows the calibration method applied to the current bracket or group. When the calibration information for the current bracket is obtained from the embedded processing method and not from a separate calibration file, the box displays *Embedded Calibration*.

For unbracketed sequences, the box displays the name of the calibration file associated with the current group in the sequence.

To change the named calibration file for an unbracketed sequence, choose File > Replace Calibration. This option is not available for bracketed sequences.

Results Grid

Figure 7 shows the results grid and its shortcut menu.

Figure 7. Results grid table and shortcut menu

| | File Name | Sample Type | Sample Name | Integration Type |
|---|------------|-------------|-------------|------------------|
| 1 | steroids02 | Standard | | Method Settings |
| 2 | steroids03 | Standard | | Method Settings |
| 3 | steroids04 | Standard | | Method Settings |
| 4 | steroids05 | Standard | | Method Settings |
| 5 | steroids13 | Standard | | Method Settings |
| 6 | steroids14 | Standard | | Method Settings |
| 7 | steroids15 | Standard | | Method Settings |
| 8 | steroids16 | Standard | | Method Settings |

Columns...
 Delete Selected Samples
 Add Sample...
 Copy Row
 Set Sorting Order...
 Send to Qual Browser

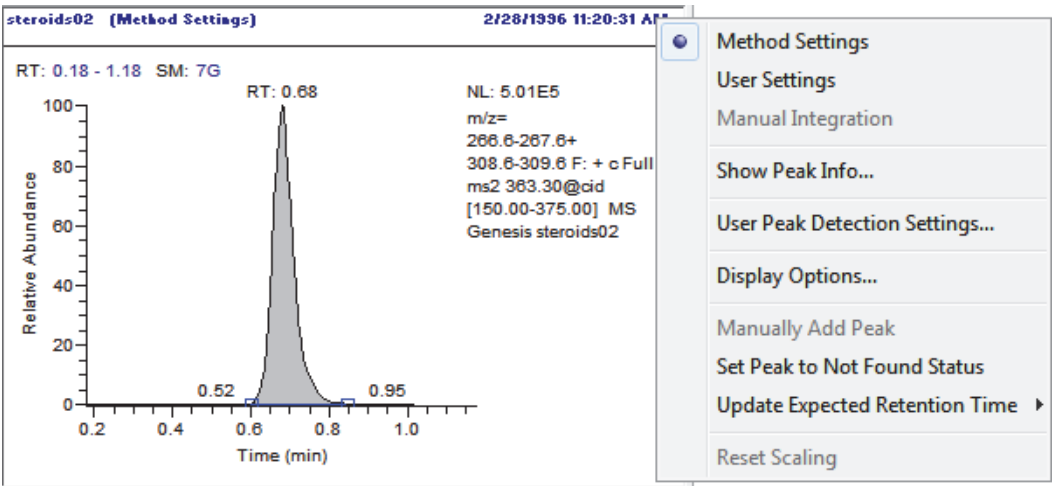
For information about the columns in this view and the shortcut menu, see “[Quan Browser Results Grid](#)” on page 84.

Chromatogram Plot and Companion Views

The chromatogram plot view displays the chromatogram for the currently selected component from the currently selected result file.

Figure 8 shows the chromatogram view and shortcut menu.

Figure 8. Chromatogram view and shortcut menu



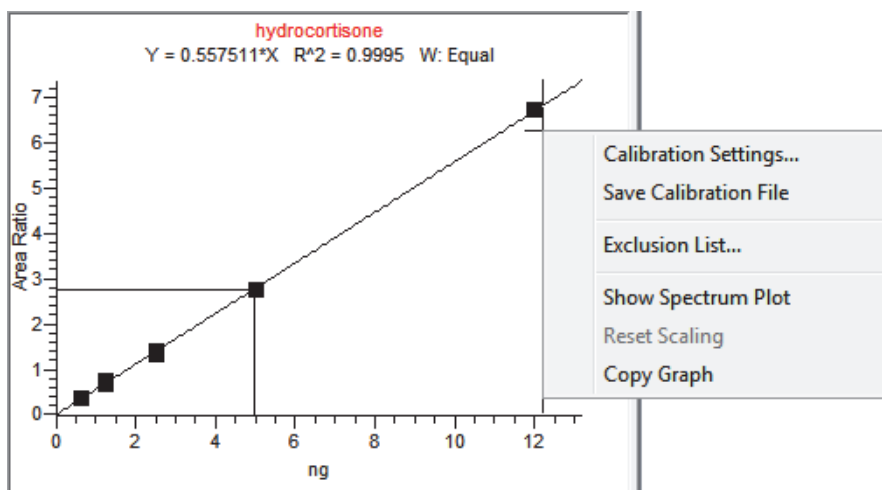
When a filter is stored within the embedded processing method for the current compound, the application applies it to the chromatogram. Adjust the chromatogram plot using the Zoom menu commands or icons in the toolbar.

The type of integration used appears in the results grid, but can be overridden. The three types are Method Settings, User Integration, and Manual Integration. Change the Integration method by using the User Identification Settings dialog box or by manually integrating a peak. For more information, see [“Using the User Identification Settings Dialog Box”](#) on page 37 and [“Integrating Chromatogram Peaks Manually”](#) on page 46.

The companion view is located to the right of the chromatogram plot view (its companion) in the lower right corner of the Quan Browser window. In the companion view you can choose either the spectrum plot view or the calibration curve view.

Figure 9 shows the calibration curve view and shortcut menu.

Figure 9. Calibration curve view and shortcut menu



Saving Changes Made in Quan Browser

In the Quan Browser window, you can do the following:

- Modify result files (RST).
- Modify processing methods (PMD).
- Create Xcalibur QuanBrowser files (XQN).
- Export the quantitation results in the results grid to a Microsoft™ Excel™ spreadsheet.
- Set up new sample sets in the results grid by adding or deleting samples; however, you cannot export the modified sample set as a new sequence file (SLD). To continue working with a modified sample set, you must save your changes to an Xcalibur Quan file (XQN).
- Evaluate manual integration settings for component peaks; however, you cannot save the results of manual integration to a result file. To continue working with manually integrated peaks, you must save your changes to an Xcalibur Quan file (XQN).

For information about changing some of the information associated with the result file, such as the sample type, sample name, or calibration level, see [“Making Changes in the Results Grid”](#) on [page 27](#). For information about changing the peak identification, detection, and integration settings, see [“Viewing and Changing Peak Identification, Detection, and Integration”](#) on [page 33](#). For information about changing the calibration settings, see [Chapter 5, “Working with the Calibration Settings.”](#)

❖ To save changes to the result files

After you make changes to result files, save the changes by choosing **File > Save All**.

❖ To save the changes to an Xcalibur QuanBrowser file

1. Choose **File > Save**.

The Save As dialog box opens.

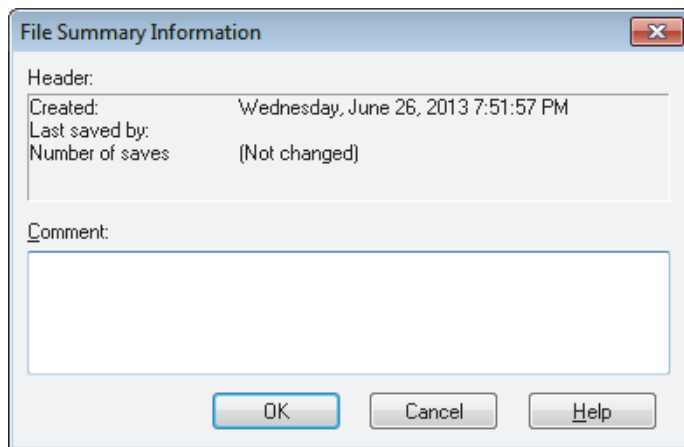
2. Select a directory in the Save In box, and type a name in the File Name box.
3. To save the file and close the Save As box, click **Save**.

❖ **To save the changes to an Xcalibur QuanBrowser file and set up an audit trail**

1. After making changes, save a file with the changes by choosing **File > Save As**.

The File Summary Information dialog box opens (Figure 10).

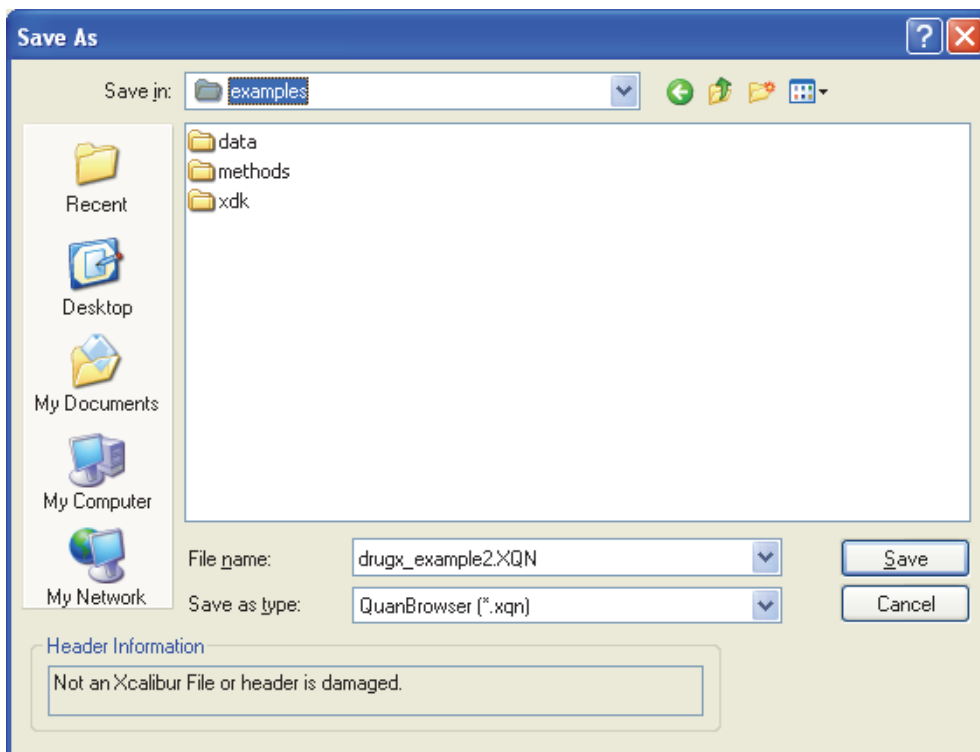
Figure 10. File Summary Information dialog box



2. Enter a comment clearly identifying the changes you made and click **OK**.

The Save As dialog box opens. The QuanBrowser (XQN) file type is the only type listed in the Save as Type list (Figure 11).

Figure 11. Save As dialog box



3. Type a name for the Xcalibur QuanBrowser file in the File Name box.
4. To save the file and close the Save As box, click **Save**.

❖ **To save user integration settings to a processing method**

1. Apply the user integration settings to one sample or to all of the samples.
2. If you applied the user integration settings for a component peak to a specific sample, and you want to save those settings to a processing method, select the sample in the results grid. If you applied the user integration to all of the samples, you do not need to select a specific sample in the results grid.
3. Choose **File > Export Method**.
The Save As dialog box opens.
4. Browse to the directory where you want to store the processing method.
5. If the selected directory folder contains processing methods, the box under the File Name box displays the method names.
6. Select the processing method that you want to modify or type a new name in the File Name box.
7. Click **OK**.

❖ **To export the information in the results grid to an Excel spreadsheet**

1. Do one of the following:
 - Choose **File > Export Data to Excel > Export Short Excel Report**.

—or—

- Choose **File > Export Data to Excel > Export Long Excel Report**.

The Excel application opens with the report displayed behind the Quan Browser window, and Quan Browser automatically saves the spreadsheet to the current working directory.

2. Save the spreadsheet.

Working with the Quan Browser Results Grid

This chapter describes how to modify the Quan Browser sequence by changing the sample type for each row, changing the level for standards and QC samples, adding and deleting rows, and changing the sort order. This chapter also describes how to use the Add Sample dialog box to change several parameter settings in a result file, including the Sample ID, Sample Name, ISTD Corr Amt, and Dilution Factor.

For information about excluding standards from the calibration curve for a component, see “Excluding Data Points from the Calibration Curve” on page 59.

Contents

- [Reviewing and Reworking the Results Grid](#)
- [Reviewing the Peak Status Column](#)
- [Making Changes in the Results Grid](#)
- [Adding and Removing Samples](#)
- [Hiding or Displaying Columns](#)
- [Changing the Sort Order](#)

Reviewing and Reworking the Results Grid

The results grid displays information about each data file in the quantitation sequence. You can review the effect of changing some of the grid parameters, including the sample type, calibration and quality control levels, internal standard correction amount, and dilution factor.

❖ To review and edit a sequence

1. To open a sequence in the Quan Browser window, do the following:
 - a. Choose **File > Open**.

The Open dialog box appears.

- b. In the File of Type list, select one of the following file types:

- Sequence List Files (SLD)
- Result Files (RST)
- Quan Browser Files (XQN)

- c. Browse to and select the file of interest and click **Open**.

When you open a bracketed sequence, the Bracket in Use list is active and displays the available brackets. The results grid displays the samples in bracket 1.

When you open an unbracketed sequence, the Group in Use list is active and displays groups broken up logically. For more information about brackets and groups, see [“Bracket/Group In Use List”](#) on [page 96](#).

When you open a result file, the application creates a one-row sequence.

2. Inspect the sequence. If the sequence contains more than one bracket, inspect each bracket. Verify that the correct samples are listed in the results grid. Ensure that each sample in the sequence is properly associated with its sample type and that the appropriate levels are associated with the standards and QC samples. Check the status in the Peak Status column as described in [“Reviewing the Peak Status Column”](#) on [page 27](#).
3. Make changes as appropriate.
- To change the sample type or the level for a standard or QC sample, follow the instructions in [“Making Changes in the Results Grid”](#) on [page 27](#).
 - To add or remove samples, follow the instructions in [“Adding and Removing Samples”](#) on [page 28](#).
 - To hide or show columns, follow the instructions in [“Hiding or Displaying Columns”](#) on [page 31](#).
 - To change the sort order of the sequence rows, follow the instructions in [“Changing the Sort Order”](#) on [page 32](#).
4. To save your changes, do the following:
- To save your edits to an Xcalibur Quan file (XQN), choose **File > Save**.

The resulting Xcalibur Quan file contains all the necessary information required to recreate the current Quan Browser session.

- To save your edits to an Xcalibur Quan file and add information to the audit trail, choose **File > Save As**.
- To update and save the result files, choose **File > Save All**.

The application updates the result files with the new information and current time stamp.

Reviewing the Peak Status Column

If the sequence contains QC samples, you can check the stability of the chromatographic method by viewing the results in the Peak Status column of the results grid in the Quan Browser window. If the calculated amount for the QC sample differs by more than the specified percentage in the processing method, the peak status is listed as QC Failed.

Note For information about setting up the QC information in the processing method and the QC levels in the sequence table, refer to the Xcalibur Data Acquisition and Processing Guide

❖ To check the stability of a sequence run that contains QC samples

1. Open the sequence of acquired and processed data in the Quan Browser window.
2. Click the **QCs** tab.
3. In the component list, select a target component.
4. View the results in the Peak Status column of the results grid.

The three possible results for a QC sample are as follows:

- **Low:** if the %Difference is < 0
- **High:** if the %Difference is > 0
- **Fail:** if the %Difference is > the user-specified percentage test value

Making Changes in the Results Grid

In the results grid, you can change the selections in the Sample Type, Level, and Integration Type columns, you can exclude or include calibration data points, and you change the text in the Sample Name column by using the Add Sample dialog box.

❖ To change the sample type

Click the **Sample Type** column to display the sample type list and select a sample type from the list.

❖ To change the level of a standard or QC sample

Click the **Level** column to display the Level list and select a sample type from the list.

❖ To exclude data points from the calibration curve

Select the check boxes in the Exclude column for each calibration standard that you want to exclude.

❖ To change the text in the Sample Name column for a sample

1. Delete the sample from the results grid by right-clicking the sample and selecting **Delete Selected Samples** from the shortcut menu.
2. Add the sample back to the results grid by following the instructions in [“Adding and Removing Samples”](#) on page 28.

❖ To change the integration type for a sample component

Do one of the following:

- To change the type from Method Settings to Manual, manually integrate a peak as described in [“Integrating Chromatogram Peaks Manually”](#) on page 46. Then, click the **Integration Type** column to display the integration type list and select **Manual** from the list.
- To change the type from Method Settings to User Integration, set up new integration settings by using the User Identification Settings dialog box. Then, click the **Integration Type** column to display the integration type list and select **User Settings** from the list.

When you select User Integration, Quan Browser integrates the component peak by using the settings in the User Identification Settings dialog box. For information about using the User Identification Settings dialog box, see [“Using the User Identification Settings Dialog Box”](#) on page 37.

—or—

- To change the type from Manual or User Integration to Method Settings, Then, click the **Integration Type** column to display the integration type list and select **Manual** from the list.

When Method Settings is selected in the Integration Type list, Quan Browser integrates the component peak by using the settings in the current processing method.

Adding and Removing Samples

To change the sample name for a sample in the results grid, delete the sample from the grid, and then add it back to the grid by using the Add Sample dialog box. For more information, see [“Add Sample Dialog Box”](#) on page 95.

❖ To add samples to the results grid

1. To minimize data entry, select a sample in the results grid that is similar to the sample or samples that you are adding.
2. Right-click the results grid to display the shortcut menu and choose **Add Sample**.

The Open Rawfile dialog box opens.

3. Browse to the appropriate folder and select the raw file or raw files to add to the sequence and click **Open**.

The Add Sample dialog box opens with the parameter settings from the data file that you selected in the results grid. By default, the Use Sample Name and Comment from Selected RawFile(s) check box is selected (Figure 12).

Tip You cannot edit the following parameters in the results grid: Sample ID, ISTD Corr Amt, Dilution Factor, Sample Name, and Comment. To enter different values for each sample, add one sample at a time.

Note After you add samples to the results grid, you can change only the Sample Type, the Level for a QC or Standard sample type, and whether a calibration standard is included or excluded from the calibration curve for a component.

Figure 12. Add Sample dialog box

The screenshot shows the 'Add Sample' dialog box with the following fields and values:

- Sample Type:** Standard (dropdown)
- Level:** 1 PG STD (dropdown)
- Sample ID:** 5 (text box)
- ISTD Corr Amt:** 0.000 (text box)
- Dilution Factor:** 1.000 (text box)
- Use Sample Name and Comment from selected rawfile(s):** ☒ (checkbox)
- Sample Name:** 1 PG STD (text box)
- Comment:** 2.5 pg/ 10 mg Internal Standard (text box)

At the bottom are buttons for OK, Cancel, and Help. A blue bracket on the right side of the dialog box groups the fields from 'Sample Type' down to 'Comment', with a line pointing to the text 'Settings from the sample selected in the results grid'.

4. Review the text entries and the Sample Type and Level selections.

5. For the Sample Name and Comments parameters, do one of the following:

- To use the information that is stored in the raw files for these parameters, keep the Use Sample Name and Comment from Selected RawFile(s) check box selected.

When you add the sample or samples to the results grid, the data system populates the Sample Name and Comments fields for the added samples with the information embedded in the raw files.

—or—

- To change the Sample Names and Comments settings for the sample or samples that you are adding, clear the Use Sample Name and Comment from Selected RawFile(s) check box. Then type the appropriate text in the Sample Name and Comments boxes.

When you add the sample or samples to the results grid, the data system populates the Sample Name and Comments fields with your text entries. When you add multiple samples at a time, the data system uses the same Sample Name and Comment for all of the added samples.

6. For the remaining parameters in the Add Samples dialog box, do the following:

- In the Sample Type list, select the appropriate sample type.

Note If the processing method associated with the selected sample in the results grid does not specify any QC levels, you cannot select the QC sample type.

- In the Levels list, select the appropriate level for standard or QC sample types.
- In the Sample ID box, type an appropriate sample ID.
- In the ISTD Corr Amt box, type a value.

The acceptable range is from 0 to 10 000.

- In the Dilution Factor box, type a value.

The acceptable range is from 0.001 to 1000.

7. Click **OK**.

The new sample or samples appear in the results grid according to the sort order specified in the Quantitation results Sorting Order dialog box. When you add a standard, Quan Browser recalculates the calibration curve for the bracket and the calculated amounts for the samples within the bracket. For information about changing the sort order, see [“Changing the Sort Order” on page 32](#) and [“Quantitation Results Sorting Order Dialog Box” on page 119](#).

8. When you finish your edits, save your changes as described in [step 4 of “Reviewing and Reworking the Results Grid” on page 25](#).

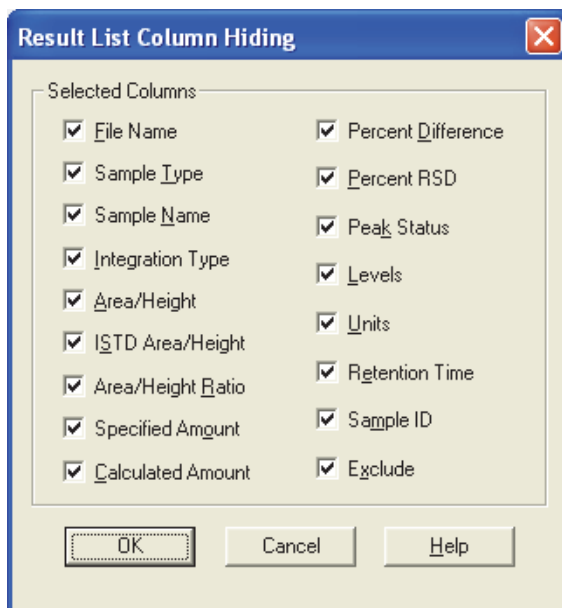
❖ **To remove samples from the sequence**

1. Select the row or rows in the sequence to delete.
2. Right-click the sequence to display the shortcut menu and choose **Delete Selected Samples**.
3. When you finish your edits, save your changes as described in [step 4](#) of “[Reviewing and Reworking the Results Grid](#)” on [page 25](#).

Hiding or Displaying Columns

The Quan Browser results grid contains up to 17 columns. Use the Result List Column Hiding dialog box ([Figure 13](#)) to display some or all of these columns.

Figure 13. Result List Column Hiding dialog box



❖ **To change the columns that the results grid displays**

1. Right-click the results grid and choose **Columns** from the shortcut menu.
The Result List Column Hiding dialog box opens.
2. Select the check box for a column heading to display the column in the results grid. Clear the check box to hide the column in the results grid.
3. Click **OK** to save your changes and close the dialog box.

Changing the Sort Order

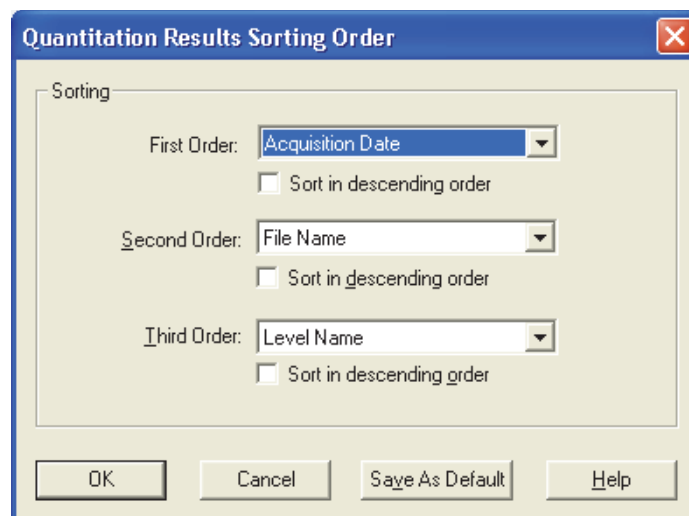
Use the Quantitation Results Sorting Order dialog box to change the order of the result data files in the results grid.

❖ **To change the sort order for entries in the results grid**

1. Right-click the results grid and choose **Set Sorting Order** from the shortcut menu.

The Quantitation Results Sorting Order dialog box opens (Figure 14).

Figure 14. Quantitation Results Sorting Order dialog box



2. To choose a heading for the primary sort of the results grid, select any of the following column headings or file properties:

- | | | |
|--------------------|---------------------|---------------|
| • <none> | • Area/Height | • Level Name |
| • %Difference | • Area/Height Ratio | • Peak Status |
| • %RSD | • Exclude | • Sample ID |
| • Acquisition Date | • File Name | • Sample Type |
| | • Integration Type | |

By default, the data system sets the first order sort to the acquisition date of the file. Select and sort with any of these sort options even if the corresponding column is not currently displayed. For example, you can sort by Sample Type, even if you have selected the Sample Name check box in the Result List Column Hiding dialog box.

3. Set any remaining column headings or file properties as the second and third sort criteria, even if the column is currently hidden.
4. To replace the default sorting criteria with your new selections, click **Save As Default**.
5. To sort the results grid, click **OK**.

Working with Peak Identification and Detection

The chromatogram plot view displays the chromatogram for the currently selected component from the currently selected result file. Most of the commands for manipulating the chromatogram plot view are available from a shortcut menu. For information about the shortcut menu, see [“Quan Browser Chromatogram Plot View”](#) on page 88.

Contents

- [Viewing and Changing Peak Identification, Detection, and Integration](#)
- [Setting the Display Options for the Chromatogram and Spectrum Plot Views](#)
- [Viewing Peak Information](#)
- [Using the User Identification Settings Dialog Box](#)
- [Integrating Chromatogram Peaks Manually](#)
- [Using the Spectrum Plot Companion View](#)
- [Reviewing Spectrum Search Results for GC/MS Data](#)

Viewing and Changing Peak Identification, Detection, and Integration

After you acquire and process data files in the Sequence Setup view, use the Quan Browser window to review the results of peak identification, detection, and integration.

❖ To review and rework peak identification, detection, and integration in Quan Browser

1. In Quan Browser, open a file (result file, sequence file, or Xcalibur Quan file), select a component in the component list, and select a sample in the results grid.
2. Right-click the chromatogram view and choose **Show Peak Info** from the shortcut menu.

The Peak Information dialog box opens.

3. Review the chromatogram peak data in these areas:

- The properties of the detected peak on the Info page
- The integration information and flags on the Flags page
- The System Suitability test results on the Suitability page
- The spectrum for the peak apex scan on the Spectrum page

For more information, see [“Viewing Peak Information”](#) on page 35.

4. Adjust the peak detection and integration in the chromatogram plot view:

- Change the detection or integration settings. See [“Using the User Identification Settings Dialog Box”](#) on page 37.
- Manually integrate peaks. See [“Integrating Chromatogram Peaks Manually”](#) on page 46.
- Change chromatogram peak labeling. To change the labels, right-click the chromatogram plot view and choose **Display Options** from the shortcut menu. In the Display Options dialog box, click the **Labels** tab. On the Labels page, select the labels to display.

5. To view spectra across the chromatogram, do the following:

- a. Display the spectrum plot view by doing one of the following:
 - From the Quan Browser menu, choose **View > Set Companion View > Show Spectrum Plot**.
 - Right-click the companion view and choose **Show Spectrum Plot** from the shortcut menu.
- b. Pin the spectrum plot view. Then, click points of interest in the chromatogram view to view the corresponding spectrum.

For more information, see [“Using the Spectrum Plot Companion View”](#) on page 48.

6. To do a detailed qualitative analysis of the chromatogram, right-click the data file of interest in the results grid and choose **Send to Qual Browser** from the shortcut menu.

The Qual Browser window opens and displays the selected result file. For more information about using Qual Browser, refer to the *Xcalibur Qualitative Analysis User Guide*.

Setting the Display Options for the Chromatogram and Spectrum Plot Views

With the exception of the Composition options for the spectrum view, the display options for the chromatogram plot and spectrum plot views in Quan Browser are the same as those for the chromatogram and spectrum views in Qual Browser.

For information about changing the display options for the chromatogram and spectrum views, refer to the *Xcalibur Qualitative Analysis User Guide*.

Viewing Peak Information

Quan Browser displays information about the currently displayed component peak, qualifier ion, or spectrum candidate in the Peak Information dialog box. The title bar contains the component name.

For more information about the Peak Information dialog box, see “[Peak Information Dialog Box](#)” on [page 108](#).

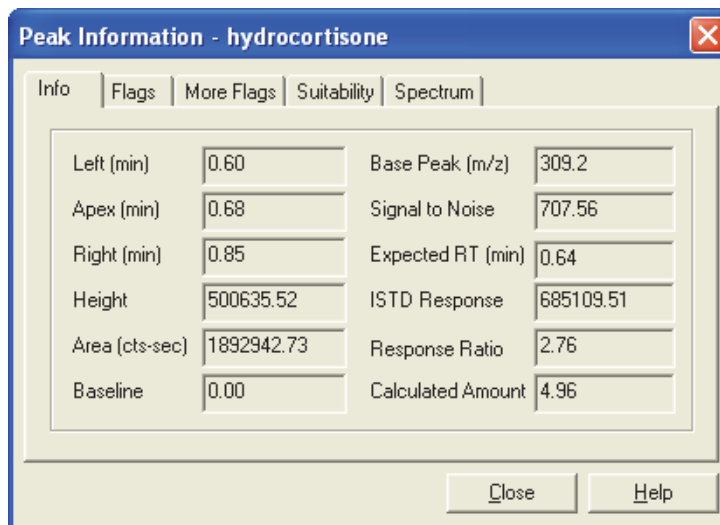
❖ To view the peak information

1. In the component list, select the component of interest.
2. Right-click the chromatogram view and choose **Show Peak Info** from the shortcut menu.

The Peak Information dialog box opens ([Figure 15](#)).

The parameters in this dialog box are read-only. If you select other components or samples, the dialog box updates the peak information for the displayed component chromatogram peak.

Figure 15. Info page – Peak Information dialog box



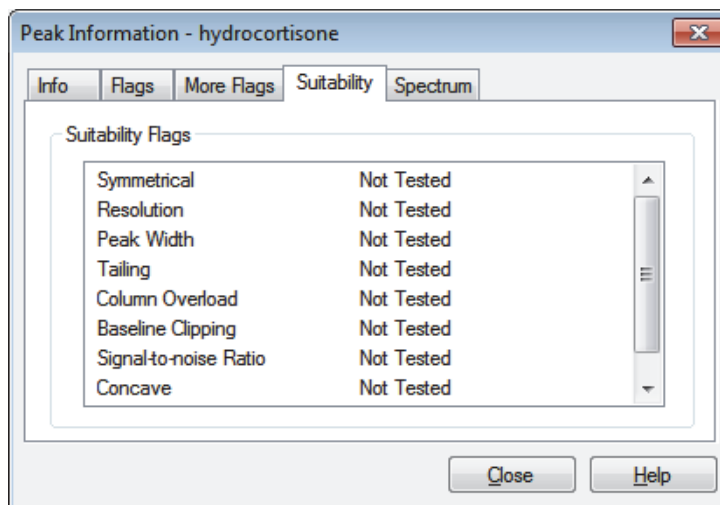
| Peak Information - hydrocortisone | | | |
|-----------------------------------|------------|-------------------|-----------|
| Left (min) | 0.60 | Base Peak (m/z) | 309.2 |
| Apex (min) | 0.68 | Signal to Noise | 707.56 |
| Right (min) | 0.85 | Expected RT (min) | 0.64 |
| Height | 500635.52 | ISTD Response | 685109.51 |
| Area (cts-sec) | 1892942.73 | Response Ratio | 2.76 |
| Baseline | 0.00 | Calculated Amount | 4.96 |

❖ To check the validity of a chromatographic peak

1. Open the Peak Information dialog box for the component of interest as described in “[To view the peak information](#)” on [page 35](#).
2. Click the **Suitability** tab.

The Suitability page opens ([Figure 16](#)).

Figure 16. Suitability page of the Peak Information dialog box



Use the parameters on the Suitability page to determine if the LC column is degrading and to identify suspicious peaks eluting at the same time as the target compound. Suspicious peaks due to highly retained compounds from a previous injection tend to have a broader than expected peak profile. Tailing peaks frequently indicate a degrading LC column.

There are three possible results for each test: Passed, Failed, or Not Tested.

To change the system suitability criteria for a component peak, you must change the criteria on the System Suitability page of the Processing Setup – Quan view, and then reprocess the raw data file or the sequence file from Sequence Setup view. For more information about setting the system suitability parameters in the processing method for a quantitative analysis, refer to the *Xcalibur Data Acquisition and Processing User Guide*.

Using the User Identification Settings Dialog Box

When you first open a sequence in Quan Browser, the application gets peak detection, calibration, and quantification information from the result file.

Within Quan Browser, use the User Identification Settings dialog box to test the effect of different parameter settings. This dialog box duplicates the parameters available on the Identification and Detection pages in the Quan view of Processing Setup.

For more information about this dialog box, see “[User Identification Settings Dialog Box](#)” on [page 129](#).

❖ To test and change the identification, detection, and integration criteria for a component peak

1. Review the displayed data for the selected component to determine if the results are consistent with your expectations:
 - Are there peaks that were not found?
 - Are neighboring peaks resolved?
 - Are tailing peaks detected properly?
2. To modify the identification and detection criteria, right-click the chromatogram plot view and choose **User Peak Detection Settings** from the shortcut menu.

The User Identification Settings dialog box opens with the Identification page displayed (see [Figure 17](#) on [page 39](#)).

3. Make changes as appropriate:
 - To change the chromatogram information or the peak detection algorithm, or to adjust the retention time window, change the settings on the Identification page.
For more information, see “[Setting the Identification Values](#)” on [page 38](#).
 - To change the detection settings, click the **Detection** tab and change the settings on the Detection page.
For more information, see “[Setting the Detection Parameters](#)” on [page 40](#).
 - If you have identified problems with noise in the peak, unresolved peaks, or peak tailing, click the **Algorithm Integration** tab and change the settings on the Integration page.

For more information about the parameters on the Integration page, see “[Integration Page – User Identification Settings Dialog Box](#)” on [page 139](#).

4 Working with Peak Identification and Detection

Using the User Identification Settings Dialog Box

- If baseline noise is interfering with peak identification or integration, modify the settings on the *Algorithm* (Genesis or ICIS) Advanced page. Use advanced options only if the standard options do not provide sufficiently selective detection criteria.

For more information about the advanced integration parameters, see [“Advanced Page – User Identification Settings Dialog Box”](#) on page 145.

- To view or change the settings on the Flags page, click the **Flags** tab.

For more information, see [“Setting the Flags Values”](#) on page 45.

4. Click **Apply** to apply the new setting to the selected sample component or **Apply to All** to apply the new settings to all of the samples in the current bracket.
5. Click **OK** to close the dialog box.
6. To save the new settings, do the following:
 - To export your settings as a new processing method, choose **File > Export Method**.
 - To save the settings in a Quan Browser file (XQN), choose **File > Save** or **File > Save As**.

Setting the Identification Values

The application uses the Identification page parameters for the following:

- Generating a chromatogram from raw data
- Identifying the component peaks in the chromatogram

For more information about this page, see [“Identification Page – User Identification Settings Dialog Box”](#) on page 130.

❖ **To set identification parameters**

1. In the User Identification Settings dialog box, click the **Identification** tab.

The Identification page opens (Figure 17).

Figure 17. Identification page – User Identification Settings dialog box

The screenshot shows the 'User Identification Settings' dialog box with the 'Identification' tab selected. The 'Name' field contains 'hydrocortisone'. The 'Plot Type' is set to 'Mass Range'. The 'Scan Filter' is '+ c Full ms2 363.30 [150.00-375.00]'. The 'Mass (m/z)' field shows '267.1, 309.1'. The 'Detector Type' is 'MS' and the 'Peak Detection Algorithm' is 'Genesis'. Under 'Retention Time', 'Expected (min)' is '0.64' and 'Window (sec)' is '30.00'. There are checkboxes for 'Use as RT Reference' and 'Adjust Using', and a 'View Width (min)' of '1.00'. At the bottom are buttons for 'OK', 'Cancel', 'Apply', 'Apply To All', and 'Help'.

2. Select the type of trace and optional trace math operation stored in the processing method in the adjacent Plot Type lists. Only certain combinations of trace types are possible.

For information about valid trace types, see “[Identification Page – User Identification Settings Dialog Box](#)” on page 130.

3. To apply a different scan filter, select a new filter from the Scan Filter list, select a new filter from the Scan Filter list and edit it, or type a new scan filter command string in the box using the scan filter format.

For information about scan filter formats, refer to the *Xcalibur Qual Browser User Guide*.

4. View or change the masses stored in the processing method.

This display area changes to accommodate the type of data required. When a single mass range is required, a single edit box displays the current value. If two mass ranges are required (as in the case of a trace defined as a Mass Range \pm Mass Range or Base Peak \pm Mass Range), this box is replaced by two boxes (in the case of Base Peak \pm Mass Range, this box is replaced by the BP and MR boxes). In the case of a TIC (no trace operator in use), analog, or digital traces, this box is blank.

4 Working with Peak Identification and Detection

Using the User Identification Settings Dialog Box

5. In the Keys box, view comments stored in the processing method. This is a read-only field.
6. Set the Retention Time parameters as follows:
 - To change the minimum width that a peak is expected to have if valley detection is enabled, type a new width in the Expected box.
 - To change the time window or to enter a new time window, type the number of seconds in the Window box.
 - To change the current view width, type the desired time in the View Width box.
7. Select a detector type from the Type list of detectors. (You configured this list using the Instrument Configuration dialog box.)
8. Select an algorithm from the Peak Detection Algorithm list and click **OK**.

The data system recalculates the current data using the specified algorithm. The application changes the default parameters for peak detection to parameters specific to that algorithm.

9. Do one of the following:
 - Click **Apply** to apply the changes to the selected sample and component in the results grid.

—or—

 - Click **Apply to All** to apply the current peak detection parameters to all samples that are currently displayed in the results grid of the Quan Browser window.

Setting the Detection Parameters

The application uses the parameters on the Detection page (see [Figure 18](#)) to confirm the identity of the component within the retention time window defined by the Identification settings. The options available on this page depend on whether the data is GC/MS data or LC/MS data.

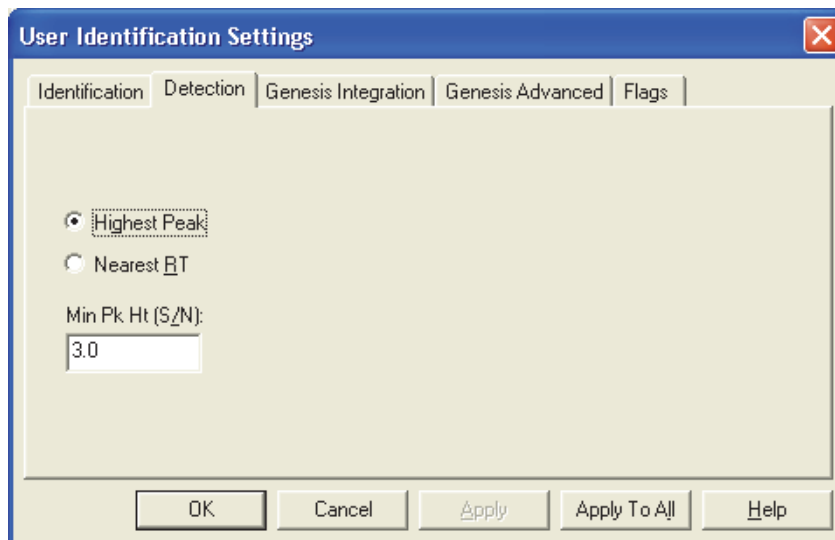
For more information about this page, see “[Detection Page – User Identification Settings](#)” on [page 134](#).

❖ **To set the detection parameters**

1. In the User Identification Settings dialog box, click the **Detection** tab.

Figure 18 shows the Detection page for LC/MS data processed with the ICIS or Genesis algorithm.

Figure 18. Detection page – User Identification Settings dialog box



2. Make changes as appropriate.

Note For information about the parameters on the Detection page for GC/MS data or for the Avalon algorithm, see “[Detection Page – User Identification Settings](#)” on [page 134](#).

3. Do one of the following:
 - Click **Apply** to apply the changes to the selected sample and component in the results grid.–or–
 - Click **Apply to All** to apply the current peak detection parameters to all samples that are currently displayed in the results grid of the Quan Browser window.

Setting the Integration Parameters

The Xcalibur data system applies the settings on the Integration page (see [Figure 19](#)) during peak integration.

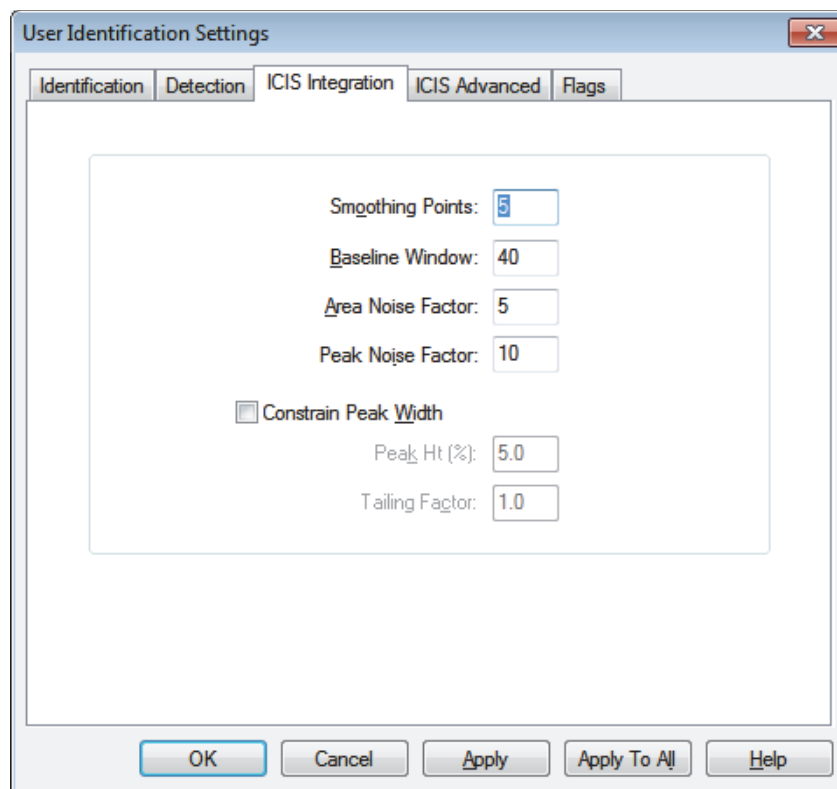
For more information about this page, see “[Integration Page – User Identification Settings Dialog Box](#)” on [page 139](#).

❖ To set the integration parameters

1. From the User Identification Settings dialog box, click the **Algorithm Integration** tab.

Figure 19 shows the ICIS Integration page.

Figure 19. ICIS Integration page – User Identification Settings dialog box



2. Make changes as appropriate.

Note For information about the parameters on the Detection page for the ICIS, Avalon, and Genesis algorithms, see “[Detection Page – User Identification Settings](#)” on [page 134](#).

3. Do one of the following:
 - Click **Apply** to apply the changes to the selected sample and component in the results grid.
 - or–
 - Click **Apply to All** to apply the current peak detection parameters to all samples that are currently displayed in the results grid of the Quan Browser window.

Setting the Advanced Parameters for the Genesis and ICIS Algorithms

The Xcalibur data system applies the Advanced page parameters (see [Figure 20](#)) during peak detection and integration.

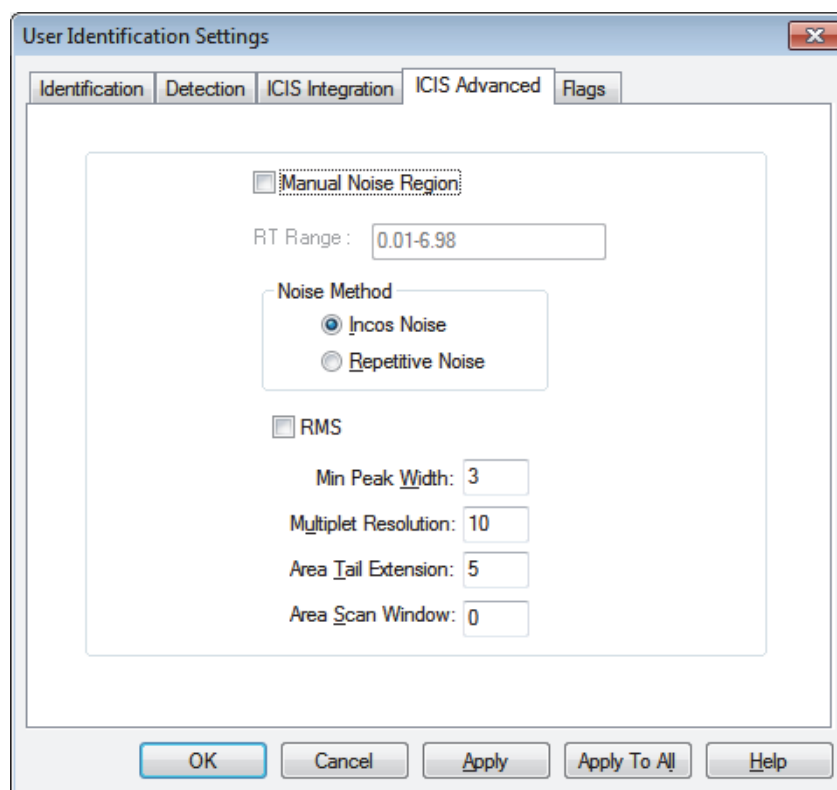
For information about the parameters on the Genesis Advanced page or the ICIS advanced page, see “[Advanced Page – User Identification Settings Dialog Box](#)” on [page 145](#).

❖ To set the advanced parameters

1. From the User Identification Settings dialog box, click the **Genesis Advanced** or the **ICIS Advanced** tab.

[Figure 20](#) shows the parameters on the ICIS Advanced page.

Figure 20. ICIS Advanced page – User Identification Settings dialog box



2. To modify the settings on the ICIS Advanced page, do the following:

a. Do one of the following:

- Select the **Manual Noise Region** check box. The Noise Method area and the RMS check box become unavailable. When you select the Manual Noise Region check box, the data system automatically calculates the RMS noise of the selected retention time range. Go to [step 2c](#).

—or—

- Select an option in the Noise Method area. The RT Range box becomes unavailable. Then, go to [step 2b](#).


b. For the Incos or Repetitive Noise options, select the **RMS** check box to use the RMS noise calculation or clear the **RMS** check box to use the peak-to-peak noise calculation.

Go to [step 2d](#).

c. To enter a retention time range for the Manual Noise Region, do one of the following:

- Type a retention time (RT) in the RT Range box.

—or—

- Click the **Manual Noise Region** icon, , in the toolbar and drag the cursor horizontally across the region of the chromatogram that you want to select as the noise region. The data system automatically fills the RT Range box with the retention time value.

Go to [step 2d](#).

d. For all of the noise methods, do the following:

- In the Min Peak Width box, type a value from **0** to **100**. This value defines the minimum number of scans required across the peak.
- In the Multiplet Resolution box, type a value from **1** to **500**. This value defines the minimum number of scans between two peaks.
- In the Area Tail Extension box, type a value from **0** to **100**. This value defines the number of scans past the peak endpoint to use in averaging the intensity.
- In the Area Scan Window box, type a value from **0** to **100**. This value defines the number of scans on each side of the peak apex.

3. Do one of the following:

- Click **Apply** to apply the changes to the selected sample and component in the results grid.

–or–

- Click **Apply to All** to apply the current peak detection parameters to all samples that are currently displayed in the results grid of the Quan Browser window.

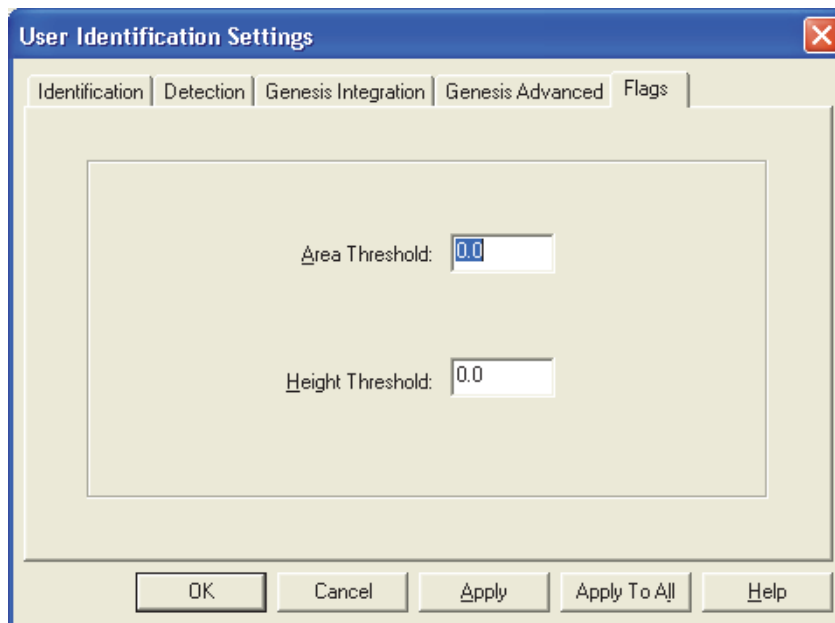
Setting the Flags Values

For more information about the Flags page, see “[Flags Page – User Identification Settings Dialog Box](#)” on [page 148](#).

❖ To use the Flags page

1. In the User Identification Settings dialog box ([Figure 21](#)), click the **Flags** tab.

Figure 21. Flags page – User Identification Settings dialog box



2. To test the validity of detected peaks, check the values on the Flags page.
3. Type a value in the current Area Threshold (AT) box. The data system sets the AT flag in the result file if the quantified peak has an area that is lower than the entered value.
4. Type a value in the current Height Threshold (HT) box. The application sets the HT flag in the result file if the quantified peak has a height that is lower than the entered value.

5. Do one of the following:

- Click **Apply** to apply the changes to the selected sample and component in the results grid.

–or–

- Click **Apply to All** to apply the current peak detection parameters to all samples that are currently displayed in the results grid of the Quan Browser window.

Integrating Chromatogram Peaks Manually

In the Quan Browser window, you can manually change the starting and ending points of chromatographic peaks.

❖ To modify and test component peak integration criteria

1. To open the sequence file that contains the calibration standards, from the Quan Browser window, choose **File > Open**.

The Open dialog box opens.

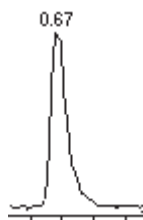
2. In the Files of Type list, select **Sequence List Files**. Browse to find the correct file. Select it and click **Open**.

The View Sample Types dialog box opens.

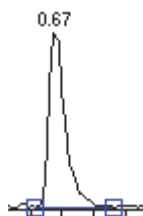
3. To accept the All Sample Types default option, click **OK**.

The data system displays the results grid, component list, chromatogram plot, and spectrum plot or calibration curve views.


4. To select a component, select the named component in the component list. The chromatogram of the component appears in the chromatogram plot view.



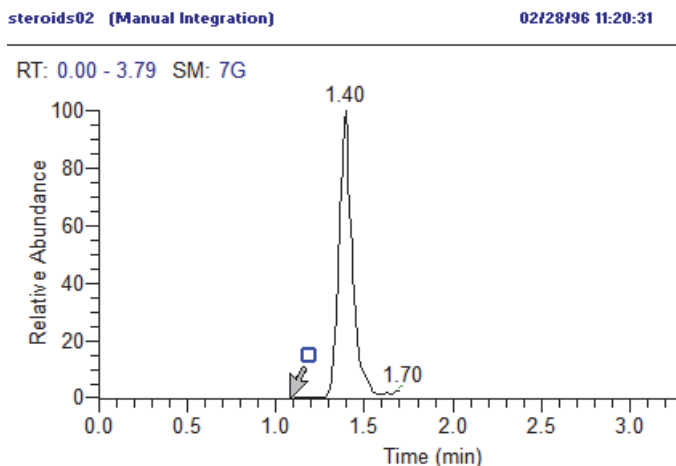
Not Found chromatogram peaks do not have blue baselines.



Found chromatogram peaks have blue baselines.

5. (Optional) Examine the chromatogram plot view as follows:
 - To replot the chromatogram with a different x axis, drag the cursor horizontally over the range that you want to expand. The application rescales the axis and replots the data with the new x -axis range.
 - To replot the chromatogram with a different y axis, drag the cursor vertically over the range that you want to expand. The application rescales the axis and replots the data with the new y -axis range.
 - To cancel the replot and return to the full range of the x axis and the highest peak normalized to 100.0 on the y axis, do one of the following:
 - Click the **Reset Scaling** icon,  in the toolbar.
 - or–
 - Choose **Zoom > Reset Scaling** from the menu bar.
6. To change the status of a peak from Found to Not Found (optional), right-click the chromatogram plot view and choose **Set Peak to Not Found Status** from the shortcut menu. The data system removes the blue baseline to the component peak, reduces the Area to 0, and changes the Peak Status to Not Found.
7. To change the status of a peak from Not Found to Found [Added] (optional), do the following:
 - a. Right-click the chromatogram plot view and choose **Manually Add Peak** from the shortcut menu.

A cursor with an arrow and a blue box appears.



- b. Drag the pointer across the desired baseline, from peak start to peak end.

The data system adds a blue baseline to the component peak, integrates the peak, displays the Area in the results grid, and changes the Integration Type to Manual Integration.

8. (Optional) Edit the baseline integration criteria of a Found or Added peak as follows:
 - a. Click the left or right square editable handle on the blue baseline of the selected peak. The data system changes the cursor to a +:



- b. Drag the handle to define a new location of the left or right peak limit. Repeat this procedure for the opposite side of the peak if required.

The data system automatically recalculates the results for the component and displays them in the results grid. If the sample is a standard, the data system replots the data and redraws the calibration curve in the calibration curve view.

Using the Spectrum Plot Companion View

The spectrum plot view displays a spectrum from the current chromatogram in the chromatogram view. View spectra from the apex, left peak edge, or right peak edge of a chromatographic peak by using commands from the shortcut menu. When the view is pinned, view scans from any part of the chromatogram by clicking the chromatogram. Initially, the data system displays the spectrum corresponding to the scan at the current chromatogram's apex retention time. If no peak is detected, the application displays the expected retention time as defined by the processing method.

Use the spectrum plot view to examine the identity of peaks and other features (such as the background) in the chromatogram.

For further analysis, including library matching of spectra, export data to Qual Browser using the Send to Qual Browser command in the shortcut menu for the results grid above the graphical views.

For more information about the spectrum plot view, see [“Quan Browser Spectrum Plot View”](#) on [page 90](#).

❖ To display the spectrum plot view

Do one of the following:

- Choose **View > Set Companion View > Show Spectrum Plot**.

—or—

- Right-click the companion view and choose **Show Spectrum Plot** from the shortcut menu.

❖ To display the spectrum for a specific time point in the current chromatogram

1. Make the Spectrum Plot view active by pinning the cell.
2. In the chromatogram view, click the time point of interest.

❖ **To use the shortcut menu commands**

Right-click the spectrum plot view and choose one of these commands from the shortcut menu:

- To display the spectrum at the apex retention time of the current chromatogram, choose **Spectrum at Peak Apex**.
- To display the spectrum at the left edge retention time of the current integration baseline, choose **Spectrum at Peak Left Edge**.
- To display the spectrum at the right edge retention time of the current integration baseline, choose **Spectrum at Peak Right Edge**.
- To change the companion view to display the calibration curve, choose **Show Calibration Curve**.
- To display the full spectrum in a normalized window, choose **Reset Scaling**.

❖ **To rescale the plot**

Use the cursor to select an area to magnify or use the Zoom menu commands and icons in the toolbar.

Reviewing Spectrum Search Results for GC/MS Data

The Xcalibur data system only performs ion spectrum searches on GC/MS data.

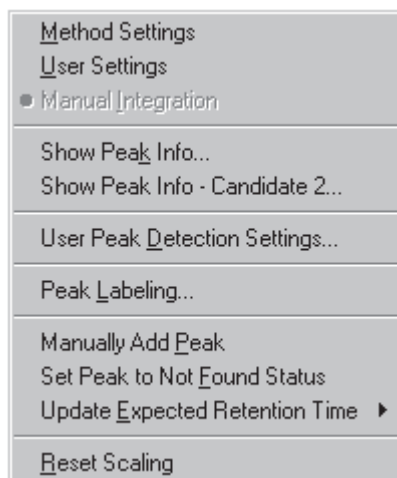
For information about setting up the spectrum search settings in the processing method, refer to the *Xcalibur Data Acquisition and Processing User Guide*.

❖ To review the results of a spectrum search in Quan Browser

1. From the Quan Browser window, choose **File > Open**. The Open dialog box opens.
2. From the Files of Type list, select either **SLD**, **RST**, or **XQN**. Browse to find the file you want to review. Select it and click **Open**. The View Sample Types dialog box opens.
3. Select the **Show All sample types** option, and click **OK**.
4. To select a component to review, click a component in the component list that has had a spectrum search as specified in the Processing Setup method. The application displays the chromatogram in the chromatogram plot view.
5. To view the plot, right-click the chromatogram plot view to display the shortcut menu.

Depending upon the spectrum search results, the data system displays 0, 1, 2, or 3 candidates in the Chromatogram Plot shortcut menu. Candidates are ranked in the order of spectrum fit as 1, 2, or 3.

This shortcut menu shows a spectrum search result that found two matching candidates.



6. Choose a spectrum search option:

- **Show Peak Info: (Not Found):** The data system did not find any peaks or could not find a spectrum match.
- **Show Peak Info:** The data system found a spectrum match and this was the best spectrum match—Candidate 1.
- **Show Peak Info - Candidate 2:** The data system found a spectrum match and this was the second best spectrum match.
- **Show Peak Info - Candidate 3:** The data system found a spectrum match and this was the third best spectrum match.

7. To review main component Candidate 1 results, choose the **Show Peak Info** command (the first of the two choices) from the Chromatogram Plot shortcut menu.

The Xcalibur system opens the Peak Information dialog box with the following title:

Peak Information – *Component*– Spectrum Candidate

and with the following pages:

Info, More Info, Flags, More Flags, Suitability, and Spectrum

Review the read-only results.

8. To review the other spectrum search results, choose the **Show Peak Info - Candidate *N*** command from the Chromatogram Plot shortcut menu.

The Xcalibur system opens the Peak Information dialog box with the following title:

Peak Information – *Component* – Spectrum Candidate

This dialog box displays the following pages:

Info, Chro, and Spectrum

The Chro page displays a total ion current plot of the Spectrum Candidate. The view is centered around the mass of interest and has the width used by the component peak display.

Review the read-only results. Repeat this step for other candidates of interest.

9. To close the Peak Information dialog box, click **Close**.

Working with the Calibration Settings

When you first open a sequence in the Quan Browser window, the Xcalibur data system performs peak detection, calibration, and quantitation according to the settings in the associated processing method.

Within the Quan Browser application, you can change the settings on the Curve, Isotope%, and Flags pages of the Calibration Settings dialog box. In addition, you can exclude one or more calibration points and one or more calibration levels from the calibration curve for each component.

Note In the Quan Browser window, you cannot change the component type (target or internal standard) or the internal standard component associated with a target compound. To change these calibration settings, you must modify the processing method in the Quan view of the Processing Setup window, and then batch reprocess the sample sequence in the Sequence Setup view. For information about creating processing methods and batch processing a sequence, refer to the *Xcalibur Data Acquisition and Processing Guide*.

Contents

- [Reviewing and Reworking the Calibration Settings in Quan Browser](#)
- [Reviewing and Reworking the Calibration Curve Settings](#)
- [Changing the Isotope Percentage Values](#)
- [Excluding Data Points from the Calibration Curve](#)

Reviewing and Reworking the Calibration Settings in Quan Browser

Use the Calibration Settings dialog box in the Quan Browser window to review and rework the calibration settings for the target components and the amounts for the internal standard components. You can use the Curve, Isotope%, and Flags pages of this dialog box to test the effect of changing the calibration curve settings, the isotope% correction values, and the threshold flags.

Use the calibration curve plot, the results grid, or the Exclusion List dialog box to exclude data points for calibration standards from the calibration curve.

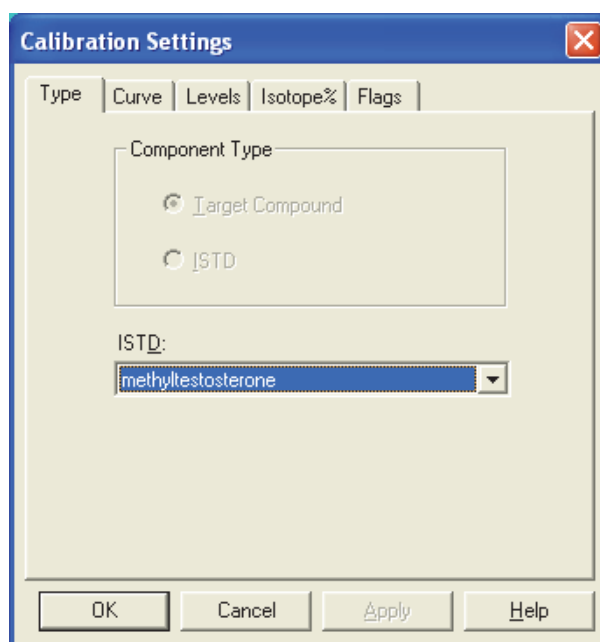
❖ To modify the calibration settings for a target component

1. Open a sequence in the Quan Browser window.
2. In the component list, select a target component.

The application automatically updates the results grid, and the chromatogram and companion views.
3. To display only calibration standards in the results grid, click the **Standards** tab.
4. If the calibration curve view is not displayed in the companion view, display it by doing one of the following:
 - Choose **View > Set Companion View > Show Calibration Curve**.
 - Right-click the chromatogram view and choose **Set Companion View > Show Calibration Curve** from the shortcut menu.
5. Inspect the calibration curve for the target component according to the criteria used in your laboratory. The calibration curve view displays the calibration equation, the goodness of fit parameter, R^2 , and the weighting, W .
6. To adjust the calibration settings, do the following:
 - a. Right-click the calibration curve view and choose **Calibration Settings** from the shortcut menu.

The Calibration Settings dialog box opens (Figure 22). For information about the pages of the Calibration Settings dialog box, see “[Calibration Settings Dialog Box](#)” on [page 98](#).

Figure 22. Calibration Settings dialog box



- b. Do the following:
 - To view the component type and the internal standard component associated with a target component, click the **Type** tab.
 - To adjust the calibration equation, weighting, or units, click the **Curve** tab and make new selections and entries on the Curve page as described in “[Reviewing and Reworking the Calibration Curve Settings](#)” on [page 55](#).
 - To view the calibration or QC levels, click the **Levels** tab. For more information, refer to the *Xcalibur Data Acquisitions and Processing User Guide*.
 - To make corrections for isotope contributions to ISTD or Target components, click the **Isotope%** tab and type new values in the boxes as described in “[Changing the Isotope Percentage Values](#)” on [page 57](#).
 - To change calibration and quantitation flag thresholds, click the **Flags** tab and type new values in the threshold boxes. For more information, refer to the *Xcalibur Data Acquisitions and Processing User Guide*.
- c. To apply any changes to the sequence, click **Apply**.
7. To exclude data points from the calibration curve, follow the instructions in “[Excluding Data Points from the Calibration Curve](#)” on [page 59](#).
8. To export the calibration settings with peak integration and detection parameters as a new processing method, choose **File > Export Method**.

Reviewing and Reworking the Calibration Curve Settings

Use the Curve page (see [Figure 23](#)) of the Calibration Settings dialog box to change the way the data system calculates and plots the calibration curve from the calibration standard data points.

For more information about the Curve page, see “[Curve Page – Calibration Settings Dialog Box](#)” on [page 98](#).

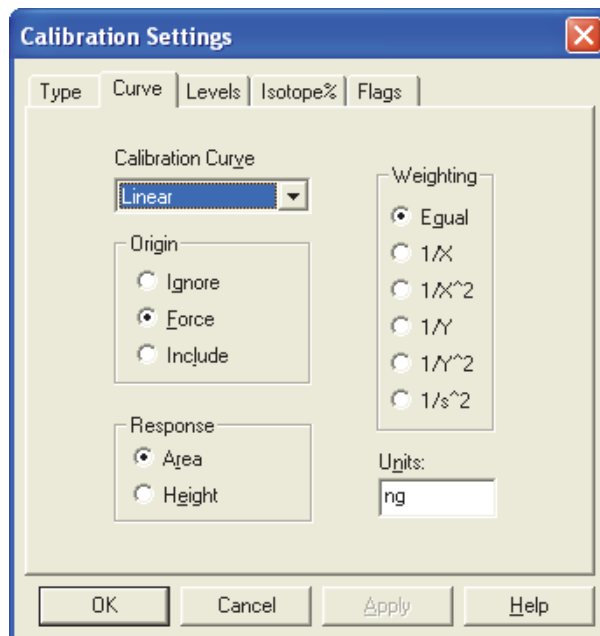
❖ To test the effect of changing the calibration curve settings for a target component

1. In the Quan Browser window, open a sequence and select the target component that you want to modify in the component list.
2. Open the Curve page of the Calibration Settings dialog box as follows:
 - a. Display the calibration curve view in the companion view by choosing **View > Set Companion View > Show Calibration Curve**.
 - b. Right-click the calibration curve and choose **Calibration Settings** from the shortcut menu.

- c. Click the **Curve** tab.

The Curve page opens (Figure 23).

Figure 23. Curve page – Calibration Settings dialog box



3. In the Calibration Curve list, select a calibration curve type:
 - If you select **Linear** or **Quadratic**, the Origin and Weighting options are available. Go to [step 4](#).
 - If you select **Point-to-Point** or **Cubic Spline**, the Weighting options are not available. Go to [step 5](#).
 - If you select **Linear Log-Log**, **Quadratic Log-Log**, **Average RF**, or **Locally Weighted**, the Origin and Weighting options are not available. Go to [step 6](#).
4. To apply weighting to the calibration points (for the Linear or Quadratic calibration curve type), select one of these options: **Equal**, **1/X**, **1/X²**, **1/Y**, **1/Y²**, or **1/s²**. This selection makes sure the data system applies the correct regression weighting method when it calculates the least-squares regression calibration curve.
5. Select how to treat the origin in the calibration curve calculation (for the Linear, Quadratic, Point-to-Point, or Cubic Spline curve type):
 - To not include the origin in the calibration curve calculation, select the **Ignore** option.
 - To require that the calibration curve passes through the origin, select the **Force** option.
 - To include the origin as one data point, select the **Include** option.

6. To select the units to be displayed on graphs and reports, type the required units label in the Units box.
7. Select how the data system determines the response:
 - To determine the response based on the integrated area of component peaks, select the **Area** option.
 - To determine the response based on the calculated height of component peaks, select the **Height** option.
8. To apply the new settings, click **Apply**.

The data system updates the values in the results grid and replots the calibration curve in the calibration curve plot view.

9. To save the new settings and close the dialog box, click **OK**.

These parameters are identical to those in the Target Compounds area on the Calibration page of Quan view in Processing Setup. These are described in more detail in “[Calibration Settings Dialog Box](#)” on [page 98](#).

Changing the Isotope Percentage Values

Use the Isotope% page (see [Figure 24](#)) of the Calibration Settings dialog box to correct data in these situations:

- An impurity in the internal standard compound that elutes at the same time as the target compound
- An impurity in the target compound that elutes at the same time as the internal standard

These parameters are identical to those in the Correction For Isotope Contribution dialog box, accessed from the Calibration page of Quan view in Processing Setup.

❖ To set the Isotope% parameters

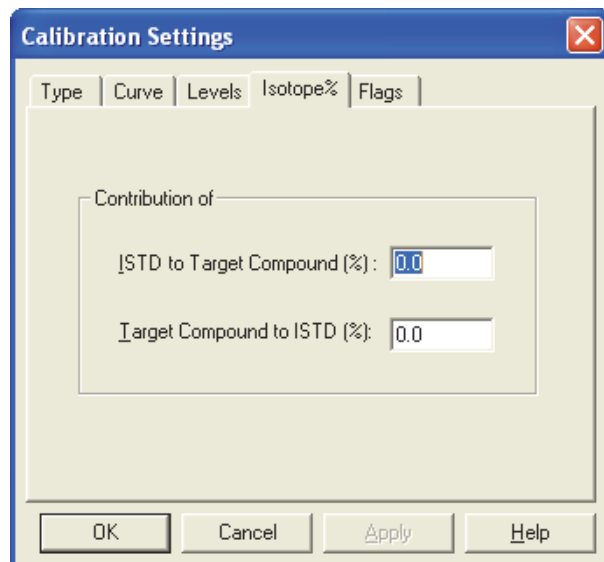
1. In the Quan Browser window, select the component of interest in the component list. Then, choose **View > Set Companion View > Show Calibration Curve** to display the calibration curve for the component.
2. Right-click the companion view and choose **Calibration Settings** from the shortcut menu.

The Calibration Settings dialog box opens.

3. Click the **Isotope%** tab.

The Isotope page opens ([Figure 24](#)).

Figure 24. Isotope% page – Calibration Settings dialog box



4. Specify the correction values as follows:

- If you have an impurity in your internal standard that elutes at the same time as the target molecule, use the ISTD to Target Compound box to type a value for the ratio $\text{ISTD}[\text{impurity}]/\text{ISTD}[\text{pure}]$, as a percentage ($\text{ratio} \times 100$).

To determine this ratio experimentally, analyze the ISTD reagent using the method to be used for quantitation of the target compound. Use the respective peak areas or heights to determine the ratio of the impurity that elutes at the same time as the target compound to the pure internal standard compound that elutes at a different retention time.

- If you have an impurity in your target compound (TM) reagent that elutes at the same time as the ISTD compound, use the Target Compound to ISTD box to type a value for the ratio $\text{TM}[\text{impurity}]/\text{TM}[\text{pure}]$, as a percentage ($\text{ratio} \times 100$).

To determine this ratio experimentally, analyze the target compound reagent using the method to be used for quantitation of the target compound. Use the respective peak areas or heights to determine the ratio of the impurity that elutes at the same time as the ISTD compound to the pure target compound that elutes at a different retention time.

The data system corrects for the impurities in the ISTD reagent, the target compound reagent, or both using the data you provide in this step, and reports the corrected amounts of ISTD and TM.

5. Click **Apply** to apply the new settings.

The data system updates the values in the results grid and replots the calibration curve.

6. To save the settings and close the dialog box, click **OK**.

Excluding Data Points from the Calibration Curve

In the Quan Browser window, you can include or exclude data points from the calibration curve in three ways:

- Click a data point in the calibration curve view and choose Include or Exclude from the shortcut menu.
- Select or clear the check boxes in the Exclude column of the results grid.
- Open the Cal Exclusion list dialog box and select the data points to be excluded or included.

Note When you include or exclude calibration standards that are shared between brackets, their status is unique to the bracket. For example, excluding a shared calibration standard in bracket 1 has no effect on its inclusion status in bracket 2.

The results grid does not display the calibration points from an external calibration file (XCAL).

IMPORTANT When you close Quan Browser, the data system erases your selections. Save your selections by choosing **File > Save As** to create a Quan Browser file (XQN).

❖ To examine the calibration curve for a target component

1. In the Quan Browser window, open the sequence of result files with the calibration curve or curves that you want to modify as follows:

- a. Choose **File > Open**.

The Open dialog box opens.

- b. In the Files of type list, select **(.sld)**. Browse to find the correct file. Select it and click **Open**.

The View Sample Types dialog box opens.


- c. Select the **Show All sample types** option and click **OK**.

The results grid, component list, chromatogram plot, and either the spectrum plot or calibration curve view open.

2. In the component list, select the component of interest.

The calibration curve for the selected component appears in the calibration curve view.

3. Examine the calibration curve data:

- To replot the data with a different x axis, drag the cursor horizontally over the range that you want to expand. The application rescales the axis and replots the data with the new x -axis range.
- To replot the data with a different y axis, drag the cursor vertically over the range that you want to expand. The application rescales the axis and replots the data with the new y -axis range.
- To cancel the replot and return to the full range for both axes, click .

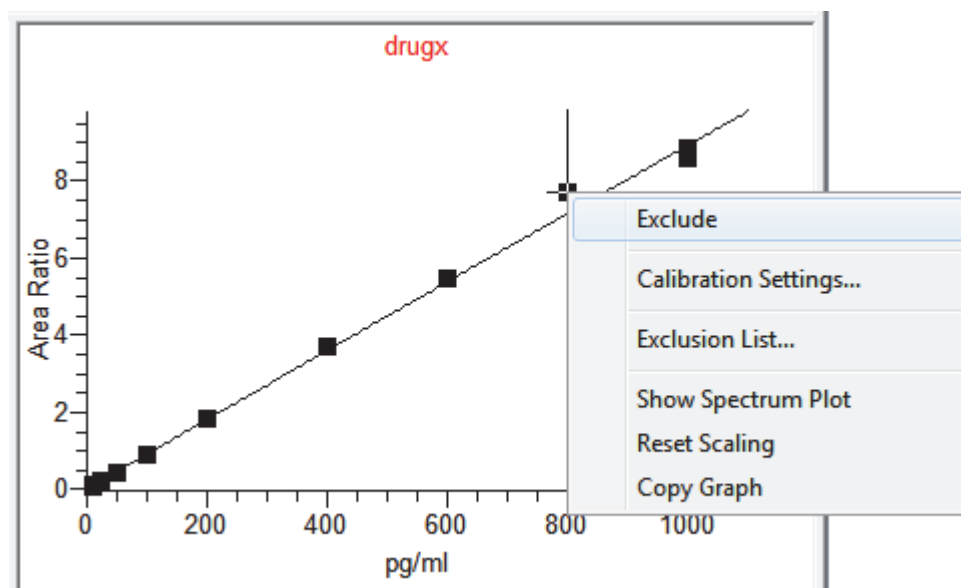
❖ **To directly include or exclude data points in the calibration curve plot view**

1. In the component list, select the target component of interest.

The calibration curve view displays the calibration curve for the selected component.

2. In the calibration curve view, right-click the point that you want to exclude and choose **Exclude** from the shortcut menu ([Figure 25](#)).

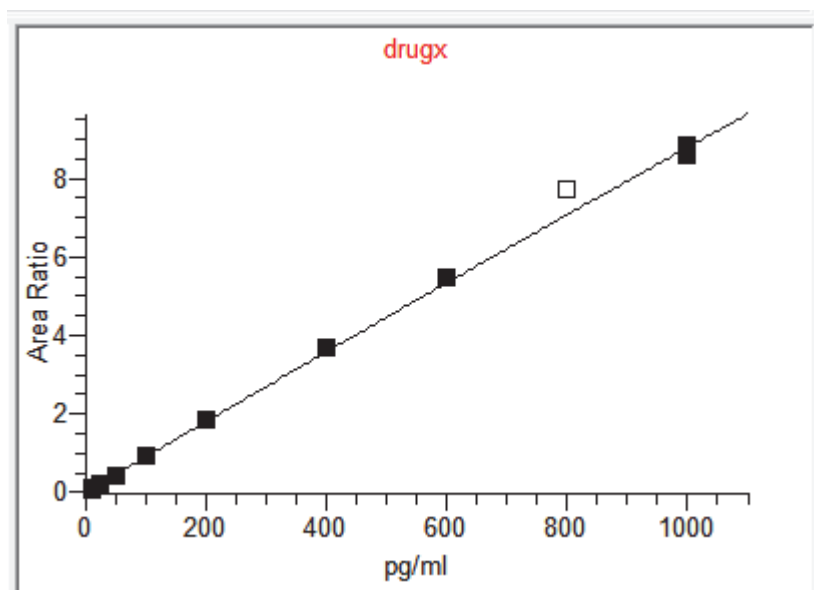
Figure 25. Selecting the Exclude command from the shortcut menu



The following occur:

- The data system recalculates the calibration curve.
- In the corresponding row of the results grid, the word “Excluded” appears in the Peak Status column and the check box is selected in the Exclude column.
- In the Cal Exclusion List dialog box, the word “Yes” appears in the Exclude column.
- In the calibration curve view, the excluded data point appears as an unfilled box ([Figure 26](#)).

Figure 26. Excluded data point appears as an unfilled box



3. To exclude an entire calibration level, exclude all of the replicate data points for the level.
4. To restore the excluded data point to the calibration curve, right-click the unfilled box and choose **Include**.

The following occur:

- The data system recalculates the calibration curve.
- In the corresponding row of the results grid, no text appears in the Peak Status column and the check box is cleared in the Exclude column.
- In the Cal Exclusion List dialog box, no text appears in the Exclude column.
- In the calibration curve view, the restored data point appears as a filled box.

❖ To include or exclude data points by using the results grid

IMPORTANT If the calibration information includes data points from an external calibration file (XCAL), the results grid does not include these data points. The results grid only includes the data points in the current sequence bracket.

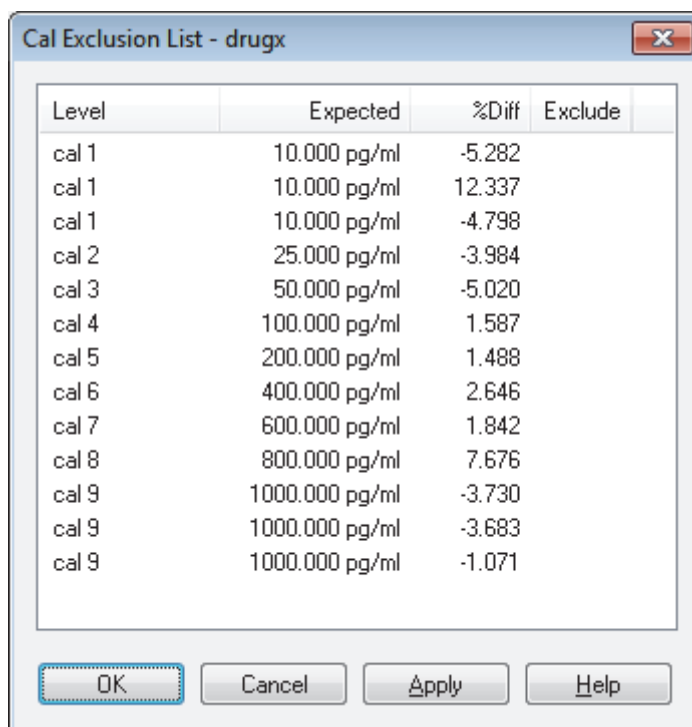
- To exclude a data point from the calibration curve, select its respective check box in the Exclude column of the results grid.
- To restore the data point to the calibration curve, clear its respective check box in the Exclude column of the results grid.

❖ **To use the Cal Exclusion List dialog box to include or exclude calibration data points**

1. In the component list, select the target component of interest.
2. Right-click the calibration curve view and choose **Exclusion List** from the shortcut menu.

The Cal Exclusion List dialog box opens (Figure 27). The Cal Exclusion List dialog box includes all the data points in the calibration curve, including the data points from an external calibration file (if used).

Figure 27. Cal Exclusion List dialog box



The dialog box lists all the replicates used in the current bracket or group and their exclusion status. The dialog box contains the following columns:

| | |
|----------|--|
| Level | Shows the name for the level. |
| Expected | Displays the expected amount for the level. |
| %Diff | Shows the percentage difference between the measured and expected amounts. |
| Exclude | Denotes excluded levels by the word <i>Yes</i> . |

3. To exclude data points and recalculate the calibration curve, do the following:

- a. Click the **Exclude** column adjacent to the data points to be excluded.
- b. Click **Apply**.

The data system does the following:

- Recalculates the calibration curve without the excluded data points.
- Updates the corresponding Peak Status and Exclude fields in the results grid to show that the data points are excluded.
- Redraws excluded data points as unfilled squares.

4. To restore excluded data points and recalculate the calibration curve, do the following:

- a. Click the **Exclude** column adjacent to the data points to be restored (on the word *Yes*).
- b. Click **Apply**.

The data system does the following:

- Incorporates the data points in the calibration and recalculates the curve.
- Updates corresponding Peak Status and Exclude fields in the results grid to show that the points are now included.
- Redraws the included data points as filled squares.

5. To close the dialog box, click **OK**.

Viewing Results and Generating Reports

Follow these procedures to review the quantitative results and generate reports.

Contents

- [Reviewing the Results of the Quantitative Analysis](#)
- [Generating Reports](#)

Reviewing the Results of the Quantitative Analysis

❖ To review and rework results before you print reports

1. Open the Quan Browser window (see [“Starting Quan Browser”](#) on page 15).

The Open dialog box opens so that you can select a file to review and rework. You can select a sequence list file (SLD), a result file (RST), or a Quan Browser file (XQN).

2. Select the file and click **OK**.

The View Sample Types dialog box opens.

3. Select either the **Show Standard and QC Sample Types** or **Show All Sample Types** option. Click **OK**.

The Quan Browser window opens with the results grid, component list, chromatogram plot view, and calibration curve or spectrum plot view.

4. In the component list, select the target component that you want to review.
5. To display calibration standards results, click the **Standards** tab.
6. Inspect the calibration curve in the calibration curve view. Evaluate the calibration curve according to the criteria used in your laboratory.
 - You can right-click the calibration curve view to display commands in the shortcut menu for modifying the information in this view.
 - You can right-click a data point in the calibration curve view to display commands in the shortcut menu for modifying the data in this view.

7. To select the first data file, click the first row in the results grid.
8. Check the entries in the results grid for peak detection and integration problems. Ensure that the selected data file corresponds to the correct level and sample type.

You can right-click the results grid to display commands in the shortcut menu for modifying the information in the results grid.

9. Change the information in any of the following columns by clicking the appropriate grid cell:
 - In the Sample Type column, click a cell and select **Standard**, **QC**, **Blank**, or **Unknown** from the list.
 - In the Integration Type column, click a cell and select **Method Settings**, **User Settings**, or **Manual Integration** from the list.
 - In the Levels column, click a cell and select another defined level from the list.
 - In the Exclude column, select or clear the **Exclude** check box in a cell to exclude or include the sample in the bracket calibration. Selecting the check box excludes the data and is indicated in the grid by *Yes*.

When a sample is shared between two brackets, you cannot change its sample type. The data system notifies you when a sample is part of two overlapping brackets if you attempt to change its Integration Type, Level, or Exclude state.

10. Inspect the component peak in the chromatogram plot view:
 - Make sure that the data system found the peak. The data system shades all found peaks gray and marks the starting and ending points with square integration markers.
 - Make sure that the data system integrated the peak properly. The shaded area should accurately represent the contribution of the component to the chromatogram.

For more information about identifying, detecting, and integrating peaks, see [Chapter 4, “Working with Peak Identification and Detection.”](#)

If necessary, perform steps [11](#) and [12](#).

11. To modify the peak detection and integration settings (optional), do the following:
 - a. Right-click the chromatogram plot view to display commands in the shortcut menu for modifying the information in this view.
 - b. Choose **User Peak Detection Settings** to display the User Identification Settings dialog box. For more information, see [“Using the User Identification Settings Dialog Box”](#) on [page 37](#).

- c. Do one or more of the following:
 - To modify the peak detection settings, click the **Detection** tab.
 - If you have problems with noise in the peak or peak tailing, click the **Integration** tab to modify the settings.
 - If baseline noise is interfering with peak identification or integration, click the **Advanced** tab to modify the settings.
12. To manually change the starting and ending points and baseline of the peak, drag the square integration markers to the desired location.
13. To review the next data file, click the next row in the results grid. Go to step 8.
14. To review and rework the QC results, do the following:
 - a. Click the **QCs** tab.
 - b. Perform steps 7 through 13 for the QCs. Evaluate the QCs according to the criteria used in your laboratory.
15. To review and rework the Blank results, do the following:
 - a. Click the **Blanks** tab.
 - b. Perform steps 7 through 13 for the blanks. Evaluate the blanks according to the criteria used in your laboratory.
16. To review and rework the Unknown results, do the following:
 - a. Click the **Unknowns** tab.
 - b. Perform steps 7 through 13 for the Unknowns. Evaluate the Unknowns according to the criteria used in your laboratory.
17. For each remaining component, select the component in the component list, and then perform steps 8 through 15.


Generating Reports

To generate reports in the Quan Browser window, follow these procedures:

- To open the Reports dialog box
- To select the samples to be included in the reports
- To set up the sample and summary reports
- To print or save the selected reports

❖ To open the Reports dialog box

In the Quan Browser window, do one of the following:

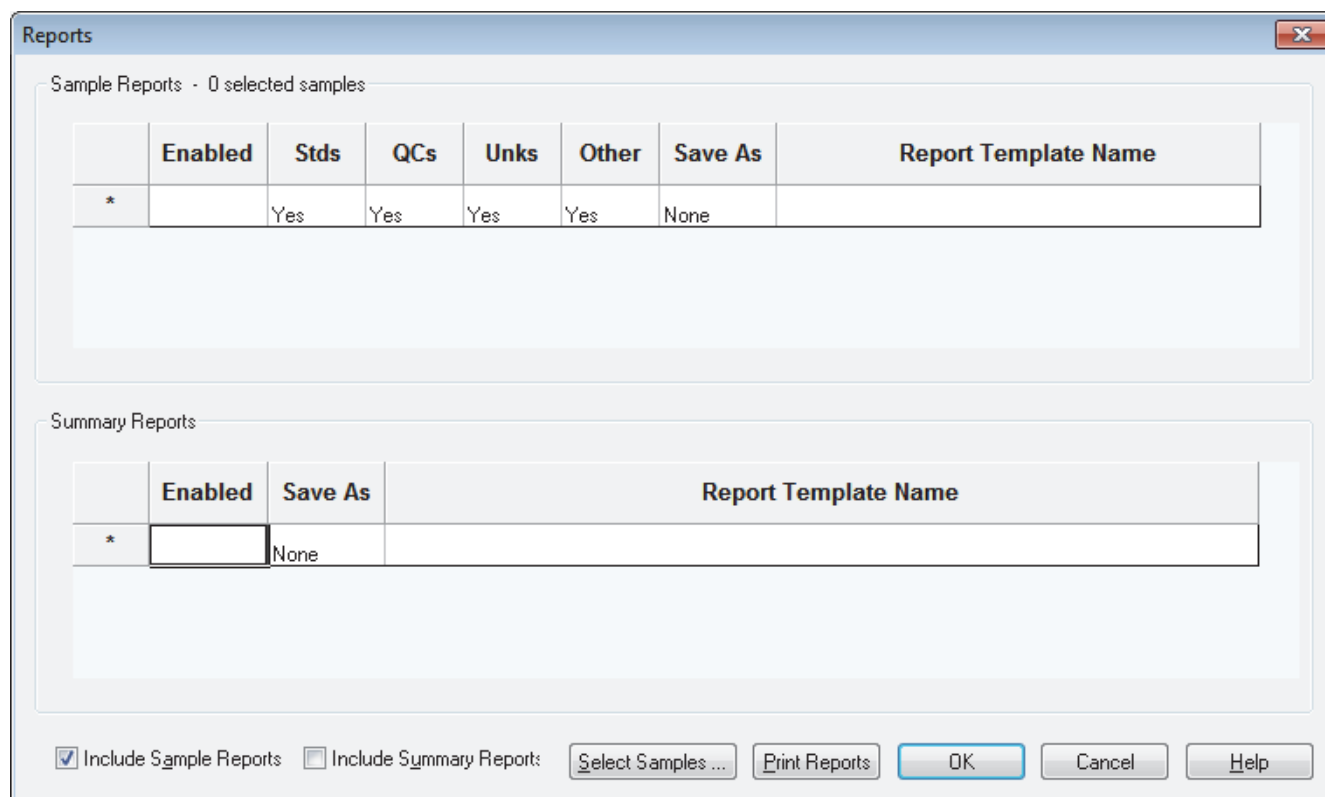
- Click the **Reports Dialog** icon, , in the toolbar.

—or—

- Choose **View > Reports Dialog** from the menu bar.

The Reports dialog box (Figure 28) duplicates the Reports view in Processing Setup. When opened, it displays the reports specified in the processing method associated with the active sequence. The displayed parameters might change as you select different brackets.

Figure 28. Reports dialog box in the Quan Browser window



Sample Reports - 0 selected samples

| | Enabled | StdS | QCs | Unks | Other | Save As | Report Template Name |
|---|---------|------|-----|------|-------|---------|----------------------|
| * | | Yes | Yes | Yes | Yes | None | |

Summary Reports

| | Enabled | Save As | Report Template Name |
|---|---------|---------|----------------------|
| * | | None | |

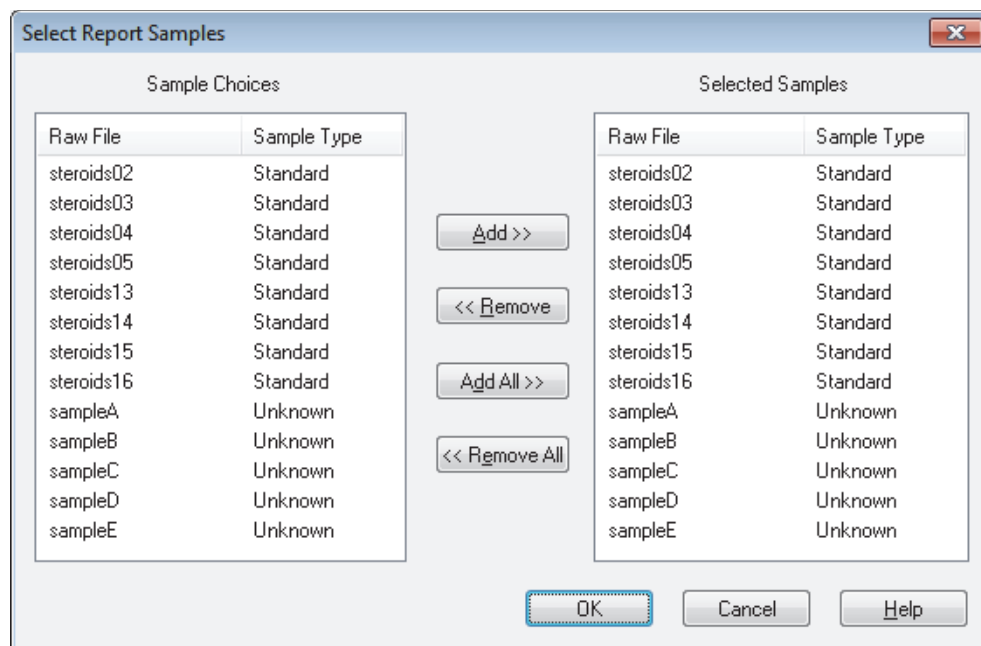
☒ Include Sample Reports
 ☐ Include Summary Reports

❖ **To select the samples to be included in the reports**

1. Click **Select Samples**.

The Select Report Samples dialog box opens. The Sample Choices box on the left lists the all of the samples in the current sequence bracket (Figure 29).

Figure 29. Select Report Samples dialog box



2. Select the samples that you want to include in the reports:

- To select a range of samples, hold the SHIFT key down. Then, click **Add** to copy the selected samples to the Selected Samples box.
- To select multiple samples, hold the CTRL key down. Then, click **Add** to copy the selected samples to the Selected Samples box.

—or—

- To include all of the samples in the reports, click **Add All**.

3. Click **OK** to close the dialog box and accept the sample list.

The Reports dialog box lists the number of samples selected.

❖ **To set up the sample and summary reports**

1. In the Sample Reports area, do the following for each sample report template that you want to use:

- a. Double-click the **Report Template Name** table cell to open the Open Template dialog box. Then, browse to the location of your report templates, select a report template, and click **Open**.

The default location of the XReport templates is as follows:

drive:\Xcalibur\templates

- b. Select the check box in the Enabled column to make the sample type columns available.

When you click another table cell, the check box is replaced with the text Yes and by default all of the sample types are selected.

- c. Click the table cell for each sample type that you want to remove from the report. and clear the check box.

The check box disappears leaving the table cell empty.

- d. In the Save As column, select one of the following;

- Select **None** if want to send the document directly to the printer.

—or—

- Select one of the available file types to save the report to the folder where your result files reside.

Figure 30 shows the selection of one sample report template for all of the selected samples. With the selection of None in the Save As column, the data system sends the reports directly to the printer.

Figure 30. Sample Reports area

| Sample Reports - 13 selected samples | | | | | | | |
|--------------------------------------|---------|------|-----|------|-------|---------|--------------------------------------|
| | Enabled | Stds | QCs | Unks | Other | Save As | Report Template Name |
| 1 | Yes | Yes | Yes | Yes | Yes | None | C:\Xcalibur\templates\QuanSimple.xrt |
| * | | Yes | Yes | Yes | Yes | None | |

2. In the Summary Reports area, do the following for each summary report that you want to use:
 - Select the check box in the Enabled column.
 - Double-click the **Report Template Name** table cell to open the Open Template dialog box. Then, browse to the location of your report templates, select a report template, and click **Open**.
3. At the bottom of the Reports dialog box, select the following:
 - To print or save the selected sample reports, select the **Include Sample Reports** check box.
 - To print or save the selected summary reports, select the **Include Summary Reports** check box.

Note If you do not select at least one of these check boxes, Quan Browser displays the following message when you click Print Reports:

No report types selected for printing.

❖ **To print or save the selected reports**

At the bottom of the Reports dialog box, click **Print Reports**.

Prior to printing reports, the data system does the following:

- Displays the Printer Activity icon at the bottom of the Quan Browser window.
- Combines the data in the sample files you selected with the report templates you selected.
- Displays the XReport Saving Report dialog box if you chose a file type (Text, Doc, HTML, PDF, RTF, or XLS) for the reports.
- Adds the report printing task to the processing queue if you chose None as the file type in the Save As column, and then displays the Printing message box.

The data system returns you to the Quan Browser window after it saves the reports to the folder where the result files reside or the reports print. The default file name format is as follows: *RawFileNameReportTemplateName*.

Quan Browser Window

Use the Quan Browser window to review sequence results for each component in each file. You can quickly review component peak identification and integration criteria. If you make any changes, you can save the new results with an audit trail describing the reason for the change.

Use Quan Browser to interactively edit processing parameters and audit changes for individual result files. It also creates new result files that contain the processing results for individual raw files. These result files include a copy of the method used to generate the results.

Result files changed using Quan Browser do not affect the original processing method. To edit processing methods, refer to the *Xcalibur Acquisition and Processing User Guide*.

Contents

- [Quan Browser Window](#)
- [Quan Browser Dialog Boxes](#)

Quan Browser Window

The Quan Browser window consists of a title bar, toolbar, menu bar, and status bar. The workspace area contains the results grid, component list, chromatogram view, and companion view. The chromatogram plot view is always displayed in the lower left portion of the window. You can display the spectrum plot view or the calibration curve view in the companion view.

For more information about the elements of the Quan Browser window, see these topics:

- “Quan Browser Title Bar” on page 75
- “Quan Browser Menu Bar” on page 76
- “Quan Browser Toolbar” on page 81
- “Quan Browser Results Grid” on page 84
- “Quan Browser Component List” on page 88
- “Quan Browser Chromatogram Plot View” on page 88
- “Quan Browser Spectrum Plot View” on page 90
- “Quan Browser Calibration Curve View” on page 92

This table provides brief descriptions of the workspace areas in the Quan Browser window.

Table 1. Quan Browser workspace (Sheet 1 of 2)

| Area | Use |
|------------------------|---|
| Results grid | View the sample rows that are retrieved from the sequence or result file. You can right-click this view to display a shortcut menu. Use this menu to review and edit the results grid. |
| Component list | View a list of all of the components contained in the current bracket. This list is in the order of user data entry in the processing method. Found/Not Found symbols are placed to the left of the component names depending on whether the Xcalibur data system found the component based on the current detection criteria for the component. |
| Chromatogram plot view | View the currently selected component chromatogram from the currently selected result file. This view, located in the lower left corner of the Quan Browser window, displays a plot of relative abundance versus time in minutes for the selected component. You can right-click this view to display a shortcut menu to review and change peak detection and integration settings. |

Table 1. Quan Browser workspace (Sheet 2 of 2)

| Area | Use |
|------------------------|--|
| Companion View | <p>Located to the right of the chromatogram plot view (its companion) in the lower right corner of the Quan Browser window. In the companion view you can choose either the spectrum plot view or the calibration curve view.</p> |
| Spectrum plot view | <p>View a plot of relative abundance (y axis) versus mass-to-charge ratio for the selected component. You can right-click this view to display a shortcut menu to display the spectrum at the left edge, apex, and right edge of the chromatogram peak. Also, you can reset the scaling of the view.</p> |
| Calibration curve view | <p>View a plot of area or area ratio (y axis) versus concentration or amount (x axis) for the selected component. You can right-click this view to display a shortcut menu to review and change calibration settings, reset the scaling of the view, and copy the graph. If you right-click a calibration point, the Calibration Point shortcut menu appears so that you can exclude/include the point, review calibration settings, reset the scaling of the view, or copy the graph.</p> |

Quan Browser Title Bar

The title bar, located in a horizontal band at the extreme top of the window, contains the following:

- Application name
- Current process
- Current file name
- Any optional parameters that are relevant

For example, if the active result file is ABC.sld, the data system might display:

Quan Browser - Browser - ABC.sld [Bracket 1, View All].

Quan Browser Menu Bar

The Quan Browser menu bar provides access to these menus:

- [File Menu – Quan Browser](#)
- [View Menu – Quan Browser](#)
- [Zoom Menu – Quan Browser](#)
- [Options Menu – Quan Browser](#)
- [GoTo Menu – Quan Browser](#)
- [Help Menu – Quan Browser](#)

File Menu – Quan Browser

This table describes the File menu commands for the Quan Browser window.

Table 2. File menu commands (Sheet 1 of 2)

| Command | Description |
|----------|--|
| Open | Open new data files. The supported file types are sequence (SLD), result (RST), and Quan Browser (XQN) files. |
| Save | Create a new Xcalibur Quan Browser file with an (.xqn) extension. This file contains all the necessary information required to recreate the current browser session. |
| Save As | Create a new Xcalibur Quan Browser file with an (.xqn) extension. This file contains all the necessary information required to recreate the current browser session. |
| Save All | Update all result files with the current information. Each result file comes from a results row in the results grid. Since each row can use method, user, or manual integration and the result file can contain only one method, the currently selected method for that row is used when creating the result file. The embedded processing method is flagged as modified. This means that each result file can potentially contain different embedded processing methods. If read back into Quan Browser, the first sample's embedded processing method is used for the method settings of the entire bracket. When read in, each sample that has a modified processing method is set in the User integration mode, and it is up to the operator to reintegrate and quantitate using a common method if desired. |

Table 2. File menu commands (Sheet 2 of 2)

| Command | Description |
|------------------------|--|
| Export Method | Export the processing method of the currently selected row. If no processing method is available, the data system uses the original processing method. |
| Export Data To Excel | |
| Short Report | Create a short Excel report. The short report displays the data contained in the current bracket and contains the same information that is available in each results grid row. |
| Long Report | Create a long Excel report. The data system reports on the data contained in the current bracket and contains the same information that is available in each results grid row, as well as other information that is not displayed in the grid. |
| Summary Information | View summary information for the current XQN file. |
| Change Dataset Name | Select a dataset from a predefined list of names. The text of this menu item might be different if the administrator chooses to use another name for a dataset. For example, this menu item might be <i>Change Job Name</i> . |
| Audit Trail | View all auditable events and changes made to data files in the current application. |
| Print Setup | Select a printer, paper, and page orientation. |
| Print | |
| Reports Dialog | Select and turn Report Templates on and off along with other options for sample reports and summary reports. |
| All Enabled Reports | Print all active and currently selected sample reports as well as all active summary reports. |
| Enable Sample Reports | Print active and currently selected sample reports. |
| Enable Summary Reports | Print all active summary reports. |
| Recently Used Files | View the paths and names of the most recently used files, located above the Exit command. The data system displays both open and closed files. Click a displayed file to load it. If the selected file is closed, the application reopens it. |
| Exit | Close the Quan Browser window. This option does not close any other Xcalibur windows. The Home Page window is not closed. |

View Menu – Quan Browser

This table describes the View menu commands for the Quan Browser window.









Table 3. View menu commands

| Command | Description |
|---------------------------|--|
| Set Companion View | |
| Show Calibration Curve | View the calibration curve view in the lower right corner of the Quan Browser window. |
| Show Spectrum Plot | View the spectrum plot view in the lower right corner of the Quan Browser window. |
| Reports Dialog | Specify the report template name to be used for sample reports or summary reports. |
| Toolbar | View or hide the toolbar. The toolbar appears if it was previously hidden or hides if it is currently displayed. |
| Status Bar | View or hide the status bar. The status bar appears if it was previously hidden or hides if it is currently displayed. |
| Show Large Toolbar | View either the large or small Quan Browser toolbar. The large toolbar is displayed when a check mark appears to the left of the command. The small toolbar is displayed when the check mark to the left of the Show Large Toolbar command is cleared. |
| Customize Toolbar | Drag any toolbar icon from this dialog box to any location on the Quan Browser toolbar and from the toolbar to this dialog box. |

Zoom Menu – Quan Browser

This table describes the Zoom menu commands for the Quan Browser window.

Table 4. Zoom menu commands

| Command | Description |
|---|--|
|  Zoom In Y | To show more detail, zoom in on the <i>y</i> axis by a factor of two (2) from the current baseline. |
|  Zoom Out Y | To show more data, zoom out on the <i>y</i> axis by a factor of two (2). |
|  Auto Range | View the chromatogram, which is normalized from the minimum to the maximum signal. (This zoom feature is recommended for PDA and UV data.) |
|  Normalize | Normalize the intensity scale of the data display to a fixed range on the <i>y</i> axis, for example, from 0–25% to 0–100%. |
|  Zoom In X | To show more detail, zoom in on the <i>x</i> axis by a factor of two (2). |
|  Zoom Out X | To show more detail, zoom out on the <i>x</i> axis by a factor of two (2) from the center. |
|  Display All | View all data on the <i>x</i> axis or all text in a report. |
|  Reset Scaling | To display the maximum amount of data, reset the scaling of both the <i>x</i> and <i>y</i> axis. |

Options Menu – Quan Browser

This table describes the Options menu commands for the Quan Browser window.





Table 5. Options menu commands

| Command | Description |
|-----------------------------|---|
| Delete <i>ComponentName</i> | Delete the currently selected component from analysis. This command results in a recalibration. |
| Masses | View or change the default settings for mass tolerance and mass precision. |
| View Stds And QCs/ View All | View either all components or just standards and quality control samples in the results grid. |
| Enable Warnings | View, as needed, all warnings boxes, even though their display has been previously suppressed. |

GoTo Menu – Quan Browser

This table describes the Go To menu commands for the Quan Browser window.

Table 6. Go To menu commands

| Command | Description |
|--|-----------------------------------|
|  Instrument Setup | Open the Instrument Setup window. |
|  Processing Setup | Open the Processing Setup window. |
|  Qual Browser | Open the Qual Browser window. |
|  Library Browser | Open the Library Browser window. |
| Xcalibur Home Page | Open the Home Page window. |

Help Menu – Quan Browser

This table describes the Help menu commands for the Quan Browser window.

Table 7. Help menu commands

| Command | Description |
|--------------------|--|
| Quan Browser Help | Open Xcalibur Help and view Help for the Quan Browser window. |
| Xcalibur Help | Open Xcalibur Help. |
| Glossary | Open the glossary. |
| How To Use Help | Open Help that describes how to use the Help viewer. |
| About Quan Browser | Open the About Quan Browser dialog box. This dialog box displays the installed version number of the Quan Browser program and the Thermo Fisher Scientific copyright notice. |

Tip You can access the manuals, as PDF files, from the Home Page – Roadmap view.

In the Home Page – Roadmap view, choose **Help > Manuals** to open a shortcut menu with links to the Xcalibur manual set.

Quan Browser Toolbar

When you install the Xcalibur data system, the Quan Browser toolbar displays these buttons.



To add or remove buttons, use the Customize Toolbar dialog box. For information about using the Customize Toolbar dialog box, see the Customize Toolbars topic in the data system Help.

This table describes the buttons and icons, by category, that can appear on the Quan Browser toolbar.

Table 8. Quan Browser toolbar buttons (Sheet 1 of 4)







| Button | Description | |
|---|---------------------------|---|
| File | | |
|  | Open | Open previously saved data files. The Xcalibur data system supports these file types: sequence (SLD), result (RST), and Quan Browser (XQN) files. |
|  | Save | When a previously saved Quan Browser file is open, clicking Save saves the recent changes to the file. When you start a new Quan Browser session, clicking Save opens the Save As dialog box, where you can save the session as a new Quan Browser file with an (.xqn) extension. This file contains all the necessary information required to re-create the current browser session. |
|  | Save As | Opens the Save As dialog box, where you can specify a name and directory location for the Quan Browser file (XQN) that contains all the necessary information required to re-create the current browser session. |
|  | Save All | <p>Update all result files with the current information. Each result file comes from a results row in the results grid. Since each row can use method, user, or manual integration, and the result file can contain only one method, the Xcalibur data system uses the currently selected method for that row when creating the result file. The system flags the embedded processing method as modified. This means that each result file can potentially contain different embedded processing methods.</p> <p>When you reopen the sequence in Quan Browser, the system uses the first sample's embedded processing method for the entire bracket's method settings. If a sample has been processed with a modified processing method, the Integration Type column of the result table displays the User Integration selection. The operator must reintegrate and quantitate using a common method.</p> |
|  | Reports Dialog | Select and turn Report Templates on or off along with other options for sample reports and summary reports. |
|  | Print All Enabled Reports | Print all active and currently selected sample reports as well as all active summary reports. |

Table 8. Quan Browser toolbar buttons (Sheet 2 of 4)















| Button | | Description |
|---|-------------------------------|---|
|  | Print Enabled Sample Reports | Print active and currently selected sample reports. |
|  | Print Enabled Summary Reports | Print all active summary reports. |
| View | | |
|  | Show Calibration Curve | View the calibration curve view in the lower right corner of the Quan Browser window. |
|  | Show Spectrum Plot | View the spectrum plot view in the lower right corner of the Quan Browser window. |
|  | Reports Dialog | Select and enable previously created report templates and other options for sample reports and summary reports. |
|  | Show Large Toolbar | View either the large or small Quan Browser toolbar. The Xcalibur data system displays the large toolbar when you select the check box to the left of the command. The system displays the small toolbar when you clear the check mark to the left of the Show Large Toolbar command. |
| Zoom | | |
|  | Zoom In Y | To show more detail, zoom in on the <i>y</i> axis by a factor of two (2) from the current baseline. |
|  | Zoom Out Y | To show more data, zoom out on the <i>y</i> axis by a factor of two (2). |
|  | Auto Range | View the chromatogram, which is normalized from the minimum to the maximum signal. (This zoom feature is recommended for PDA and UV data.) |
|  | Normalize | Normalize the intensity scale of the data display to a fixed range on the Y-axis, for example, from 0–25% to 0–100%. |
|  | Zoom In X | To show more detail, zoom in on the <i>x</i> axis by a factor of two (2). |
|  | Zoom Out X | To show more detail, zoom out on the <i>x</i> axis by a factor of two (2) from the center. |
|  | Display All | View all data on the <i>x</i> axis or all text in a report. |
|  | Reset Scaling To Full Scale | To display the maximum amount of data, reset the scaling of both the <i>x</i> and <i>y</i> axis. |

Table 8. Quan Browser toolbar buttons (Sheet 3 of 4)









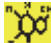


| Button | Description | |
|---|----------------------------|--|
| Noise | | |
|  | Manual Noise Region | <p>Specify a noise region manually.</p> <p>❖ To select a manual noise region</p> <p>Click the Manual Noise Region icon, , and drag the cursor horizontally across the region of interest in the chromatogram.</p> <p>The data system marks the region with a red baseline and updates noise value in the Signal to Noise box on the Info page of the Peak Information dialog.</p> <p>The application calculates noise based on the data points you select. It uses all selected data points as noise points and calculates noise based on those points. You can select the noise region from an individual trace or different noise regions from multiple traces.</p> <p>To make this button active, open a raw file and select a chromatogram.</p> |
|  | Delete Manual Noise Region | <p>Remove a designated manual noise region.</p> <p>❖ To remove a manual noise region</p> <ol style="list-style-type: none">1. Click the Delete Manual Noise Region icon, , and drag the cursor over the region that was previously selected as the noise region.2. Release the mouse button to delete the noise region. |
| Options | | |
|  | View Std's And QC's | View all components or just standards and quality control samples in the results grid. |
| GoTo | | |
|  | Instrument Setup | Open the Instrument Setup window. |
|  | Processing Setup | Open the Processing Setup window. |
|  | Qual Browser | Open the Qual Browser window. |
|  | Library Browser | Open the Library Browser window. |
|  | Xcalibur Home Page | Open the Home Page window if it is closed or display the Home Page window if it is already open. This command closes the Instrument Setup window so that all instrument setup methods are closed when samples that use these methods are run from the Home Page window. |

Table 8. Quan Browser toolbar buttons (Sheet 4 of 4)

| Button | Description |
|---|---|
| Help | |
|  Quan Browser Help | View Help for Quan Browser. Opens the Xcalibur Help system to the Quan Browser section. |

Quan Browser Results Grid

The results grid ([Figure 31](#)) displays sequence or result sample data entries. Each row defines a result file and its associated parameters. Selecting any cell within the results grid causes the additional views to refresh, using data from the newly selected sequence row. The Xcalibur data system validates any changes made to the sequence row and updates the chromatogram plot, calibration curve, and spectrum plot views automatically.

Figure 31. Results grid

| | File Name | Sample Type | Sample Name | Integration Type | Area |
|----|------------|-------------|-------------|------------------|---------|
| 1 | steroids02 | Standard | | Method Settings | 1892943 |
| 2 | steroids03 | Standard | | Method Settings | 956506 |
| 3 | steroids04 | Standard | | Method Settings | 531348 |
| 4 | steroids05 | Standard | | Method Settings | 243328 |
| 5 | sampleA | Unknown | | Method Settings | 1862850 |
| 6 | sampleB | Unknown | | Method Settings | 905012 |
| 7 | sampleC | Unknown | | Method Settings | 5453768 |
| 8 | sampleD | Unknown | | Method Settings | 278722 |
| 9 | sampleE | Unknown | | Method Settings | 489300 |
| 10 | steroids13 | Standard | | Method Settings | 1859643 |
| 11 | steroids14 | Standard | | Method Settings | 884533 |
| 12 | steroids15 | Standard | | Method Settings | 515035 |
| 13 | steroids16 | Standard | | Method Settings | 4629083 |

In the Options menu, if you choose to view to View Stds and QCs, the results grid displays three tabs: All, Stds, and QCs. The All page displays all Stds and QCs, the Stds page displays only standards, and the QCs page displays only QCs.

In the Options menu, if the viewing preference is set to View All, the results grid displays five tabs: All, Stds, QCs, Blanks, Unknowns. The All page in this case displays all Stds, QCs, Blanks and Unknowns. The Stds page displays only standards, the QCs page displays only QCs, the Blanks page displays only blanks, and the Unknowns page displays only unknowns.

Depending on the current settings of the Result List Column Hiding dialog box, the results grid can display different parameters for each sample row.

These topics describe the columns in the results grid view and the shortcut menu:

- [Results Grid Columns](#)
- [Results Grid Shortcut Menu](#)

Results Grid Columns

This table describes the columns in the results grid of the Quan Browser window.

Table 9. Results grid parameters (Sheet 1 of 2)

| Column | Description |
|--|---|
| File Name (read-only) | View the raw file that contains the acquisition data for this run. |
| Sample Type | <p>View or change the sample type that has been assigned to the sample. Every sample must be designated one of these four basic types: Standard, QC, Blank, or Unknown.</p> <p>You cannot change sample type of a sample that is shared between two brackets. If you try to make this type of change, the data system displays a warning message:</p> <p style="padding-left: 40px;">You cannot change the sample type of a sample shared between brackets. Reverting back to original sample type.</p> |
| Integration Type | <p>View or change the method the used to integrate the peak.</p> <p>The selections are Method Settings, User Integration, and Manual Integration.</p> <ul style="list-style-type: none"> • Select Method Settings to apply the integration parameters defined during the processing setup phase. • Select User Integration to override the integration parameters within Quan Browser. • Select Manual Integration when the integration baseline was selected by dragging the baseline handles. |
| Area Or Height (read-only) | View the integrated area (count-seconds) or height (counts) under the detected peak. |
| Internal Standard (ISTD) Area/Height (read-only) | View the integrated area (count-seconds) or height (counts) under the detected Internal Standard peak. |
| Area Ratio or Height Ratio (read-only) | View the area or height ratio between the selected peak and the Internal Standard peak. |
| Specified Amount (read-only) | View the amount of the component at the Cal or QC level. |
| Calculated Amount (read-only) | View the component amount, as determined by the response ratio and calibration curve. |
| % Difference (read-only) | View the percentage difference between the calculated amount and the specified amount. |

Table 9. Results grid parameters (Sheet 2 of 2)

| Column | Description |
|--|--|
| % Relative Standard Deviation (read-only) | View the standard deviation of the difference between the calculated amount and the specified amount, as expressed as a percentage of the specified value. |
| Peak Status (read-only) | View the peak status: Low: if the %Difference is < 0 High: if the %Difference is > 0 Fail: if the %Difference is > the QC fail percentage test value |
| Level | View or change the calibration or QC level of the sample. To change this value, click the grid cell and select another level from the list. |
| Units (read-only) | View the units used for quantity or concentration. |
| Retention Time (read-only) | View the retention time in minutes at the peak maximum. |
| Sample ID (read-only) | View the unique sample identification string. |
| Exclude | Indicate whether to include or exclude the sample point from the calibration curve. Include or exclude a data point by clicking the grid cell; select the check box to include the data point or clear it to exclude the data point. |

Results Grid Shortcut Menu

This table describes the menu commands in the shortcut menu for the results grid in the Quan Browser window.

Table 10. Results grid shortcut commands

| Command | Description |
|-------------------------|--|
| Column | <p>Open the Result List Column Hiding dialog box, where you can modify the columns that appear in the results grid.</p> <p>For more information, see “Result List Column Hiding Dialog Box” on page 126.</p> |
| Delete Selected Samples | Remove the currently selected samples from the results grid view. If you delete standard samples, the data system recalibrates the data. If you delete either standard samples (stds) or quality control samples (QCs), the application recalibrates the (%) Relative Standard Deviation. |
| Add Sample | Select a new file from a browse dialog box. The data system adds it to the results grid view and sorts it according to the sort order. If you add standard samples, the application recalibrates the data. If you add either standard samples (stds) or quality control samples (QCs), the Xcalibur system recalibrates the (%) Relative Standard Deviation. |
| Copy Row | Duplicate the currently selected row and add it to the next row of the results grid view. If the row added is a standard sample (std), then a recalibration takes place. If it is either a standard sample (std) or a quality control sample (QC), then the data system recalculates the (%) Relative Standard Deviation. |
| Set Sorting Order | Set the sort order for the samples in the results grid view. |
| Send To Qual Browser | Open the result file for the sample that you selected in Qual Browser. |

Quan Browser Component List

The component list on the right side of the window lists the components in the processing method associated with the currently loaded sequence. To select a component of interest, select the component name from the list. All Quan Browser views are automatically updated to reflect the settings and data for the selected component. The component list provides the only method of selecting a component of interest.

Figure 32. Component list for the drugx example set



Quan Browser Chromatogram Plot View

The chromatogram plot view is the graphical display pane in the lower left portion of the Quan Browser window.

This table describes the shortcut menu commands for the chromatogram plot view.

Table 11. Chromatogram plot view shortcut commands (Sheet 1 of 2)

| Command | Description |
|--------------------|--|
| Method Settings | Reset the integration method to those specified in the original method stored within the result file. |
| User Settings | Apply a set of user-defined peak detection parameters to the integration of this peak. See “Using the User Identification Settings Dialog Box” on page 37 for more information. |
| Manual Integration | If active, this command indicates that the integration baseline has been adjusted by dragging the control boxes on the plot. This command is inactive until you adjust the baseline by using the graphical dragging operation at least once. |
| Show Peak Info | View peak information in a read-only format. The information displayed reflects the settings for the currently displayed chromatographic peak. Select other compounds or other result files to get detailed information on the found peak. To close the dialog box, click Close . The compound identification name is displayed in the title bar. For example: Peak Information – drugx |

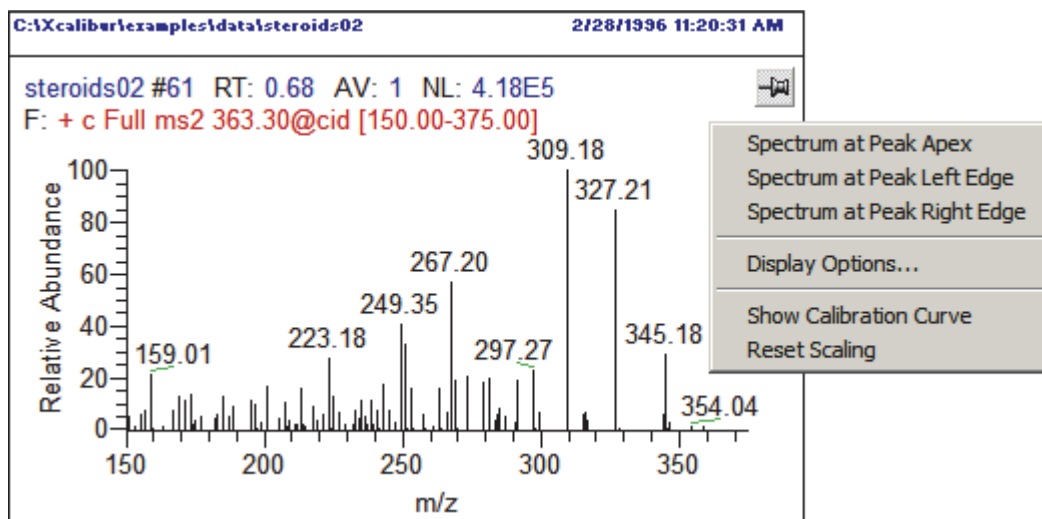
Table 11. Chromatogram plot view shortcut commands (Sheet 2 of 2)

| Command | Description |
|--------------------------------|--|
| User Peak Detection Settings | Apply unique peak detection parameters to the chromatogram. The tabbed pages on this dialog box differ for GC and LC processing methods. These differences are described in the appropriate sections. The type of processing used to create the result file comes from the processing method (also stored within the result file) and are not available for changes within Quan Browser. |
| Display Options | Place labels, such as retention time, scan number, base peak, signal-to-noise, flags, area, and height, on your chromatograms. This menu command is not active if no chromatogram data is displayed in the chromatogram plot view. |
| Manually Add Peak | Manually place a baseline on the Chromatogram Plot and set the current integration method for the current component to Manual Integration. The Xcalibur data system changes the entry in the Integration Type box of the results grid view. If no peak has been detected for the currently selected compound, there is no integration baseline on the chromatogram plot for you to manually adjust. |
| Set Peak To Not Found Status | Tag the current peak as Not Found, do another peak search (possibly with new integration parameters), or do a manual integration to restore the peak. |
| Update Expected Retention Time | Update the retention time specified in the processing method with the retention time that the Xcalibur data system detected. Update Component In Current Row: This command updates the retention time of the current component in the current row of the active sequence and recalculates the data for the row in the results grid view. Update All Components In Current Row: This command updates the retention time of all the components in the current row of the active sequence and recalculates the data for the row in the results grid view. |
| Reset Scaling | This command resets the plot scale to include the full peak in a normalized window. |

Quan Browser Spectrum Plot View

When you select the spectrum plot view (Figure 33) located in the lower right corner of the Quan Browser window as the companion plot, the spectrum plot view displays a spectrum from the peak apex of the current chromatogram.

Figure 33. Spectrum plot view with shortcut menu displayed



To display the spectrum plot view in the companion view, do one of the following:

- In the toolbar, click the **Show Spectrum Plot** icon, .

—or—

- From the menu bar, choose **Set Companion View > Show Spectrum Plot**.

You can select a spectrum from a different time point in the chromatogram in two ways:

- Pin the spectrum plot view, and then click a time point in the chromatogram.
- Right-click the spectrum plot view and select **Spectrum At Peak Left Edge** or **Spectrum At Right Edge** from the shortcut menu.

This table describes the shortcut menu commands for the spectrum plot view. For more information about working with the spectrum plot view, see [“Using the Spectrum Plot Companion View”](#) on [page 48](#).

Table 12. Spectrum plot view shortcut commands

| Command | Description |
|-----------------------------|--|
| Spectrum At Peak Apex | View the spectrum at the apex retention time of the displayed chromatogram peak. |
| Spectrum At Peak Left Edge | View the spectrum of the left edge of the displayed chromatogram peak. The left edge is defined at the retention time corresponding to the left baseline box (handle). |
| Spectrum At Peak Right Edge | View the spectrum of the right edge of the displayed chromatogram peak. The right edge is defined at the retention time corresponding to the right baseline box (handle). |
| Display Options | Select the style, color, labels, axis, and normalization options for your spectrum from the Display Options dialog box. This menu command is not active if no spectrum data is displayed in the spectrum plot view. |
| Show Calibration Curve | View the calibration curve view in the lower right corner of the Quan Browser window. |
| Reset Scaling | Reset the plot scale to include the full peak in a normalized window. |

Quan Browser Calibration Curve View

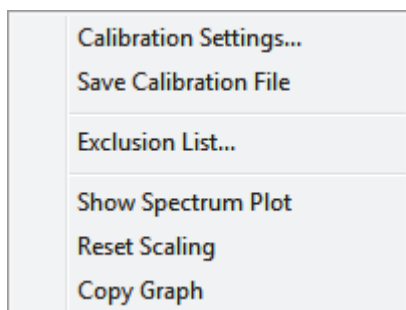
To display the calibration curve view in the companion view, do one of the following:

- In the toolbar, click the **Show Calibration Curve Plot** icon, .

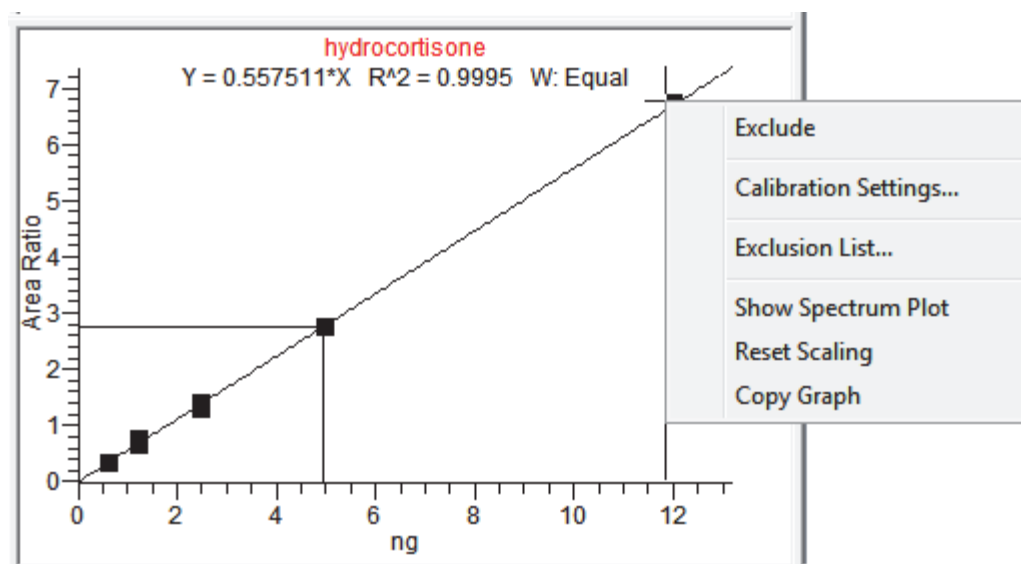
—or—

- From the menu bar, choose **View > Set Companion View > Show Calibration Curve**.

Right-click in the calibration curve view (except on a calibration point) to display the shortcut menu for the view.



Right-click a calibration point on the calibration curve in the calibration curve view to display the following shortcut menu.



This table describes the shortcut menu commands for the calibration curve view.

Table 13. Calibration curve view shortcut commands

| Command | Description |
|-----------------------|--|
| Exclude | If the data point is currently included in the calibration, selecting the Exclude command forces the calibration curve to be recalculated without the selected data point. The excluded entry in the results grid view is updated to show that this point is excluded and the excluded data point is redrawn on the calibration curve using an unfilled square. |
| Include | <p>If the current data point was previously excluded, the first menu item is Include. Select this command to include the data point and recalculate the calibration curve. The data system updates the excluded entry in the results grid to show that this point is now included and redraws the included data point on the calibration curve using a filled square.</p> <p>Note You can include or exclude samples that are shared between brackets. Sample status is unique to the bracket—that is, excluding a shared sample in bracket 1 has no effect on its inclusion status in bracket 2.</p> |
| Calibration Settings | Select values for the following parameters: type, curve, levels, isotope%, and flags. These options appear when you display a valid calibration curve. |
| Save Calibration File | Name and save the calibration file to the directory of your choice. |
| Exclusion List | Open the Exclusion List dialog box and select points in the calibration curve to include or exclude. |
| Show Spectrum Plot | View the spectrum plot view in the lower right corner of the Quan Browser window. |
| Reset Scaling | Reset the plot scale to include the full peak in a normalized window. |
| Copy Graph | Copy the calibration curve view to the clipboard so that you can transfer it to another open application, such as Microsoft Word, using the Paste command. This operation is useful when you are writing a report and want to include the calibration curve. |

Quan Browser Dialog Boxes

The Quan Browser window has the following dialog boxes:

- “Add Sample Dialog Box” on page 95
- “Bracket/Group In Use List” on page 96
- “Cal Exclusion List Dialog Box” on page 97
- “Calibration Settings Dialog Box” on page 98
- “Display Options Dialog Box in Quan Browser” on page 107
- “Masses Dialog Box” on page 107
- “Peak Information Dialog Box” on page 108
- “Quantitation Results Sorting Order Dialog Box” on page 119
- “Reports Dialog Box” on page 121
- “Result List Column Hiding Dialog Box” on page 126
- “Select Level Dialog Box” on page 127
- “Select Report Samples Dialog Box” on page 128
- “User Identification Settings Dialog Box” on page 129
- “View Sample Types Dialog Box” on page 149

Add Sample Dialog Box

Use the Add Sample dialog box to change information about the sample or samples that you are adding to the results grid in the Quan Browser window.

For information about adding samples to the results grid and changing the settings in the result files for the parameters listed below, see [“Adding and Removing Samples”](#) on [page 28](#).

This table describes the parameters in the Add Sample dialog box.

Table 14. Add Sample dialog box parameters

| Parameter | Description |
|--|---|
| Sample Type | Select a sample type for a new sample. |
| Level | Select a level for a new sample. |
| Sample ID | Enter the sample ID for a new sample. |
| Sample Name | Enter the sample name for a new sample. |
| ISTD Corr Amt | Enter the ISTD amount for a new sample. |
| Dilution Factor | Enter the dilution factor for a new sample. |
| Use Sample Name and Comment from Selected RawFile(s) | To edit the Sample Name and Comments fields, clear this check box. To use the Sample Name and Comment embedded in the raw file, select this check box. |
| Comment | Enter a comment about the new sample. |

Bracket/Group In Use List

Use this list to select the current bracket or group being browsed. Select a new bracket or group from the list to refill the results grid view with the samples from the selected bracket. The Xcalibur data system updates all additional plots and dialog boxes automatically.

This table describes the Bracket/Group In Use list that is located above the results grid of the Quan Browser window.

Table 15. Bracket/Group In Use list parameters

| Parameter | Description |
|----------------|--|
| Bracket in use | If a bracketed sequence is open, the Bracket in use list is active and displays the available brackets numbered sequentially from one to the number of brackets in the sequence, for example: Bracket 1, Bracket 2, Bracket 3, and so on. If the sequence contains multiple brackets, the data system selects the first bracket in the list and displays the samples related to that bracket in the results grid view. If the sequence contains only one bracket, then the list displays only one entry: Bracket 1. |
| Group in use | <p>If an unbracketed sequence is open, the Group in use list is active and displays groups broken up logically. For example: Group 1, Group 2, Group 3, and so on. Groups are separated by standards and are created using the following rules:</p> <p>The data system scans the sequence file in chronological order until the first standard (either Std Clear or Std Update) is encountered. The first standard begins a new group. The application continues looking through the sequence, adding all samples to this group until the next standard after the first nonstandard is found. There are two deviations from this rule. The first exception is that the first group does not have to start with a standard. The first sample, by definition, begins the first group. The second exception is that any Std Clear begins another group even if it immediately follows a Std Update.</p> |

Cal Exclusion List Dialog Box

In the Quan Browser window, you can exclude a data point from the calibration curve in three ways. You can right-click a data point in the calibration curve view and select Exclude from the shortcut menu. You can select the appropriate check box in the Exclude column of the results grid. Or, you can use the Cal Exclusion List dialog box. For more information about excluding data points from the calibration curve, see [“Excluding Data Points from the Calibration Curve”](#) on page 59.

The Cal Exclusion List dialog box displays a list of all the data points used in creating the current calibration curve. When you use an external calibration file (.xcal) with an unbracketed sequence, the results grid does not include the data points from the external calibration file. The results grid only includes the data points from the current sequence bracket. The data points from the external calibration file are accessible in the Cal Exclusion List dialog box or the calibration curve plot.

The title bar of the Cal Exclusion List dialog box contains the name of the selected component:

Cal Exclusion List – *ComponentName*

The exclusion list displays all the replicates used in the current bracket or group and their exclusion status. If there is a *Yes* in the exclude column, then the Xcalibur data system excludes this replicate data point in the calibration curve plot. If the column is blank, then the Xcalibur system includes the replicate in the calibration curve plot. To change the status of any replicate, click the exclude column in the row containing the replicate data point. This changes the value. Click **Apply** to make the changes, but keep the Exclusion List dialog box visible. Click **OK** to accept the changes and close the dialog box. Click **Cancel** to abandon the changes and close the dialog box.

This table describes the parameters in the Cal Exclusion List dialog box.

Table 16. Cal Exclusion List dialog box parameters

| Parameter | Description |
|-----------|---|
| Level | View the read-only calibration level names for the sample data points displayed in the calibration curve view. |
| Expected | View the read-only expected quantity values for the data points displayed in the calibration curve view. |
| % Diff | View the read-only percent difference values for the data points displayed in the calibration curve view. These values are the percentage difference between the calculated amount and the expected amount. |
| Exclude | Click each row (replicate) in the Exclude column to turn Yes on and off. Yes indicates that the data point is excluded from the data points displayed in the calibration curve view. |

Calibration Settings Dialog Box

The Xcalibur data system displays particular pages in the Calibration Settings dialog box, based on whether the component is an Internal Standard or a Target. For more information, see, “[Reviewing and Reworking the Calibration Settings in Quan Browser](#)” on page 53.

For an internal standard component, the following page is available:

- “[Type Page – Calibration Settings Dialog Box](#)” on page 105

For a target component, the following pages are available:

- “[Curve Page – Calibration Settings Dialog Box](#)” on page 98
- “[Flags Page – Calibration Settings Dialog Box](#)” on page 100
- “[Isotope% Page – Calibration Settings Dialog Box](#)” on page 101
- “[Levels Page – Calibration Settings Dialog Box](#)” on page 104
- “[Type Page – Calibration Settings Dialog Box](#)” on page 105

Curve Page – Calibration Settings Dialog Box

Use the Curve page to modify component type settings that you specified on the Calibration page in the Quan view of the Processing Setup window.

To test the results of the new setting, click **Apply** or **OK**.

This table describes the parameters on the Curve page of the Calibration Settings dialog box.

Table 17. Curve page parameters – Calibration Settings dialog box (Sheet 1 of 2)

| Parameter | Description |
|---|---|
| Calibration Curve | View or change the selected calibration curve type. |
| Origin | |
| Specify how to use the origin in your calibration curve. You can choose to ignore the origin as a data point, force the curve to pass through it, or include it as a single point on the calibration curve. | |
| Ignore | Exclude the origin as a valid point in your calibration curve. If you select this option, the calibration curve might or might not pass through the origin. |
| Force | Specify that the calibration curve passes through the origin of the data point plot. |
| Include | Include the origin as a single data point in the calculation of the calibration curve. If you select this option, the calibration curve might or might not pass through the origin. |

Table 17. Curve page parameters – Calibration Settings dialog box (Sheet 2 of 2)

| Parameter | Description |
|------------------|---|
| Response | |
| | Use the area or the height of the target compound peak to acquire the data used for the calibration. |
| Area | Use the area of the target compound peak to acquire the data used for the calibration. |
| Height | Use the height of the target compound peak to acquire the data used for the calibration. |
| Weighting | |
| Equal | Give all calibration data points equal weight during the least-squares regression calculation of the calibration curve. |
| 1/X | Specify a weight value of 1/X for all calibration data points during the least-squares regression calculation of the calibration curve. The data system assigns a weight value of the inverse of their quantity. |
| 1/X ² | Specify a weight value of 1/X ² for all calibration data points during the least-squares regression calculation of the calibration curve. The data system assigns a weight value of the inverse of the square of their quantity. |
| 1/Y | Specify a weight value of 1/Y for all calibration data points during the least-squares regression calculation of the calibration curve. The data system assigns a weight value of the inverse of their response (or response ratio). |
| 1/Y ² | Specify a weighting of 1/Y ² for all calibration data points during the least-squares regression calculation of the calibration curve. The data system assigns a weight value of the inverse of the square of their response (or response ratio). |
| 1/s ² | Specify a weighting of 1/s ² for all calibration data points during the least-squares regression calculation of the calibration curve. Calibrants at a given level are weighted by the inverse of the standard deviation of their responses (or response ratios). For this weighting factor to be used, there must be two or more replicates at each level. If only one calibrant is available for any level, 1/s ² weighting cannot be used. |
| Units | |
| | View or change the units set on the Calibration page of the Quan view of Processing Setup. The units are also used in reports and in Quan Browser. |

Flags Page – Calibration Settings Dialog Box

The Flags page of the Calibration Settings dialog box displays the current calibration and quantitation flags in use for the selected compound. The Xcalibur data system uses these values in determining if the calibration or quantitation is within user specified criteria. They do not alter the way calculations are made. An entered value of zero forces the flag to be false.

During entry of the values in the Quantitation Flags box, the Xcalibur data system checks to make sure that the relationships between the four fields are maintained. If an entry in one field forces a change to occur in another field, the Automatic Adjustment message box is displayed with the following message:

The last change made has forced an automatic change in one or more of the remaining edit boxes. This is due to the fact that each value and the acceptable ranges for that value are linked.

This table describes the parameters on the Flags page of the Calibration Settings dialog box.

Table 18. Flags page parameters – Calibration Settings dialog box

| Parameter | Description |
|---------------------------|--|
| Calibration Flag | |
| R-Squared | View or change the current value of R-squared. The data system sets the R-Squared (RS) flag in the result file if the computed coefficient of determination is lower than the entered R-Squared value. |
| Quantitation Flags | |
| Limit Of Detection | View or change the current value for the limit of detection. The data system sets the Limit of Detection (LOD) flag in the result file if the peak concentration is less than the entered Limit of Detection. |
| Limit Of Quantitation | View or change the current value for the limit of quantitation. The data system sets the Limit of Quantitation (LOQ) flag in the result file if the peak concentration is less than the entered Limit of Quantitation. |
| Linearity Limit | View or change the current value for the linearity limit. The data system sets the Limit of Linearity (LL) Flag in the result file if the peak concentration is greater than the entered Linearity Limit. Linearity Limit + 10 percent is the calibration curve's <i>x</i> -axis upper value (default) when displayed in the Quan Browser window. The user can then go on to change this <i>x</i> -axis range. |
| Carry Over Limit | View or change the current value for the carryover limit. The data system sets the Carry Over Limit (COL) flag in the result file if the peak concentration is greater than the entered Carry Over Limit. |

Isotope% Page – Calibration Settings Dialog Box

Use the Isotope (%) page of the Calibration Settings dialog box to correct for one or both of the following:

- An impurity in the internal standard compound that elutes at the same time as the target compound
- An impurity in the target compound that elutes at the same time as the internal standard compound

❖ To determine the percent impurity in the internal standard reagent

Analyze the ISTD reagent using the method to be used for quantitation of the target compound. Use the respective peak areas or heights to determine the ratio of impurity to pure compound.

❖ To change the percent impurity for the internal standard reagent,

Type a new value in the Contribution of ISTD to Target Compound box, as a percentage (ratio \times 100%).

The valid range is 0.00 to 100.00 percent.

❖ To determine the percent impurity in the target compound reagent

Analyze the Target reagent using the method to be used for its quantitation. Use the respective peak areas or heights to determine the ratio of impurity to pure compound.

The valid range is 0.00 to 100.00 percent.

❖ To change the percent impurity for the target compound reagent

Type a new value in the Contribution of Target Compound to Internal Standard box, as a percentage (ratio \times 100%).

The valid range is 0.00 to 100.00 percent.

This table describes the parameters on the Isotope (%) page of the Calibration Settings dialog box.

Table 19. Isotope (%) page parameters – Calibration Settings dialog box (Sheet 1 of 2)

| Parameter | Description |
|-----------------------------|--|
| Contribution of | |
| ISTD To Target Compound (%) | <p>Specifies the percent impurity in the internal standard reagent:</p> $\text{Percent impurity in ISTD reagent} = \frac{\text{ISTD [impurity]}}{\text{ISTD [pure]}} \times 100$ <p>Where:</p> <p>ISTD [impurity] is an impurity compound in the internal standard reagent that elutes at the same time as the target compound.</p> <p>ISTD [pure] is the pure internal standard compound.</p> <p>The valid range is 0.00 to 100.00 percent.</p> <p>The data system uses this percentage (converted to a ratio) as the value for x in the following impurity correction expressions:</p> $\text{ISTD [corr]} = [\text{ISTD [obs]} - y \text{ Target [obs]}] / [1 - yx]$ $\text{Target [corr]} = [\text{Target [obs]} - x \text{ ISTD [obs]}] / [1 - yx]$ <p>Where:</p> <p>ISTD [corr] is the corrected amount of the internal standard compound.</p> <p>ISTD [obs] is the apparent amount of the internal standard compound, as measured by the application at the retention time for the internal standard compound. This peak consists of ISTD [corr] + Target [impurity].</p> <p>Target [corr] is the corrected amount of the target compound.</p> <p>Target [obs] is the apparent amount of the target compound, as measured by the application at the retention time for the target compound. This amount consists of Target [corr] + ISTD [impurity].</p> <p>See the parameter description “Target Compound To ISTD” on page 103 for a complete description of variable y.</p> |

Table 19. Isotope (%) page parameters – Calibration Settings dialog box (Sheet 2 of 2)

| Parameter | Description |
|----------------------------|---|
| Contribution of | |
| Target Compound To ISTD | <p>Specifies the percent impurity in the target compound:</p> $\text{Percent impurity in Target reagent} = \frac{\text{Target [impurity]}}{\text{Target [pure]}} \times 100$ <p>Where:</p> <p>Target [impurity] is an impurity compound in the target compound reagent that elutes at the same time as the internal standard.</p> <p>Target [pure] is the pure target compound.</p> <p>The data system uses this percentage (converted to a ratio) as the value for y in the following impurity correction expressions:</p> $\text{ISTD [corr]} = [\text{ISTD [obs]} - y \text{ Target [obs]}] / [1 - yx]$ $\text{Target [corr]} = [\text{Target [obs]} - x \text{ ISTD [obs]}] / [1 - yx]$ <p>Where:</p> <p>ISTD [corr] is the corrected amount of the internal standard compound.</p> <p>ISTD [obs] is the apparent amount of the internal standard compound, as measured by the data system at the retention time for the pure internal standard compound, ISTD [pure]. This peak consists of ISTD [corr] + Target [impurity].</p> <p>Target [corr] is the corrected amount of the target compound.</p> <p>Target [obs] is the apparent amount of the target compound, as measured by the data system at the retention time for the pure target compound, Target [pure]. This amount consists of Target [corr] + ISTD [impurity].</p> <p>See the parameter description “ISTD To Target Compound (%)” on page 102 for a complete description of variable x.</p> |

Levels Page – Calibration Settings Dialog Box

Use the Levels page of the Calibration Settings dialog box to review the current component type settings that you specified in the processing method. These are read-only boxes.

Table 20. Levels page parameters – Calibration Settings dialog box

| Parameter | Description |
|---|---|
| Cal Level | View the calibration levels for the selected component. |
| Amount | View the amount of target compound used for each calibration level. |
| QC Level/ Amount/%Test Table settings | View QC (quality control) level names, quality control level amounts, and %test values. QC samples containing known amounts of a component can be utilized to help ensure the accuracy of an analysis. The data system measures the quantity of the QC component in the same manner as unknown components and compares the measured quantity with a user-defined expected quantity and a user-defined percent test. |
| QC Level | View the quality control levels for the selected component. The application can accommodate up to 15 QC levels. |
| Amount | View the quantity used for each QC level, as defined by the user. |
| % Test | View a value for the acceptable difference (as a percent) between the known amount and the calculated (measured) amount of each QC level. |
| Units | View the units set on the Calibration page of the Quan view in Processing Setup. The application also uses the units in reports and in Quan Browser. |

Type Page – Calibration Settings Dialog Box

The parameters on the Type page of the Calibration Settings dialog box changes depending on whether the selected component is an internal standard component or a target component.

For an internal standard component, use the Type page to change the Amount and Units settings that you specified on the Calibration page in the Quan view of the Processing Setup window.

For an analysis that contains more than one internal standard component, use the Type page to associate a different internal standard component with the selected target component.

This table describes the parameters for an internal standard component on the Type page of the Calibration Settings dialog box.

Table 21. Internal Standard component parameters

| Parameter | Description |
|---|--|
| Component Type | |
| The data system recognizes two component types for component calibration purposes: Target Compound components and Internal Standards (ISTD) components. You can initially define, select, and characterize components of interest using the Calibration page in the Quan view of the Processing Setup window. | |
| In the Quan Browser window, you can change the amount of the internal standard component and the display units. The options in the Component Type area are unavailable. | |
| Target Compound | For an internal standard component, this option is not selected. |
| ISTD | For an internal standard component, this option is selected. |
| Internal Standard (ISTD) | |
| Amount | Specifies the amount of the selected internal standard component that is added to each sample. To change the amount, type amounts with up to three decimals of precision in the Amount box. |
| Units | Specifies the units used for the internal standard amount or concentration, for example, ng or pg/mL. You can change the units. |

This table describes the parameters for target component on the Type page of the Calibration Settings dialog box.

Table 22. Target component parameters

| Parameter | Description |
|--|---|
| Component Type | |
| <p>The data system recognizes two component types for component calibration purposes: Target Compound components and Internal Standards (ISTD) components. Components of interest are initially defined, selected, and characterized using the Calibration page in the Quan view of the Processing Setup window. These parameters can then be revised using the Type page of the Calibration Settings dialog box available in the Quan Browser window.</p> | |
| Target Compound | For a target component, this option is selected. |
| ISTD | <p>For a target component, this option is not selected.</p> <p>The ISTD option is unavailable if you processed the data with an external standard calibration.</p> |
| ISTD | <p>This list contains all the internal standard components defined in the processing method.</p> <p>If the processing method contained more than one internal standard component, you can select a different internal standard component for the selected target component from the list.</p> |

Display Options Dialog Box in Quan Browser

In the Quan Browser window, use the pages of the Display Options dialog box to select Style, Color, Labels, Axis, and Normalization settings. The parameters on these pages depend on whether the view in the active cell is a chromatogram or a spectrum.

For information about the parameters on these pages, refer to the *Xcalibur Qualitative Analysis User Guide* or the Xcalibur Help.

Note The parameters on the Style, Color, Labels, Axis, and Normalization pages of the Dialog Options dialog box depend on whether the view in the active cell is a chromatogram or a spectrum. The same pages for the chromatogram and spectrum views are available from the real-time display view and the Qual Browser, Quan Browser, and Processing Setup windows.

Masses Dialog Box

Use this dialog box to specify tolerance and precision settings for the mass data displayed in the chromatogram, spectrum, map, and ion map plots.

Specify the default values for tolerance and precision on the Mass Options page of the Xcalibur Configuration dialog box.

This table describes the parameters in the Masses dialog box.

Table 23. Masses dialog box parameters

| Parameter | Description |
|-----------------------|--|
| Mass tolerance | |
| Mass tolerance | Specify the value for mass tolerance. Type a value in the range of 0.1 to 50 000 and select units to apply to the value. The Xcalibur data system uses the tolerance value to create the mass range limit. |
| Units | Specify the units of measurement in which the data system processes your data. Select mmu (millimass units) or ppm (parts per million). |
| Mass precision | |
| Decimals | Specify the number of decimal places (mass precision) that the data system uses to display mass values. You can specify from 0 to 5 decimal places. The number of decimal places applies to the mass data in a window. |

•

Peak Information Dialog Box

The Peak Information dialog box provides read-only information about the currently displayed chromatogram peak (for GC/MS or LC/MS data) or one of the peaks used by the spectrum search or ion ratio confirmation routines (for GC/MS data). You can use this dialog box to quickly get detailed information on the found peak by selecting other components or other result files.

The current component name is displayed in the title bar, for example:

Peak Information – *drugx*

Depending on the peak type, the Peak Information dialog box displays one or more of these pages:

- “No Peak Page – Peak Information Dialog Box” on page 109
- “Info Page – Peak Information Dialog Box” on page 110
- “Flags Page – Peak Information Dialog Box” on page 111
- “More Flags Page – Peak Information Dialog Box” on page 114
- “Chro Page – Peak Information Dialog Box” on page 115
- “Info or More Info Page – Peak Information Dialog Box” on page 116
- “Spectrum Page – Peak Information Dialog Box” on page 118
- “Suitability Page – Peak Information Dialog Box” on page 118

For more information about reviewing the information in the Peak Information dialog box, see “[Viewing Peak Information](#)” on page 35.

[Table 24](#) describes the possible peak types. [Table 25](#) lists the pages displayed in the Peak Information dialog box for each peak type.

Table 24. Peak types for the Peak Information dialog box (Sheet 1 of 2)

| Peak type | Description |
|--------------------|---|
| No peak | <p>If no peak has been found for the component, the Xcalibur data system displays only the No Peak page with the following message:</p> <p>No Peak Found. Cannot show Peak Info.</p> |
| Standard peak (LC) | <p>A standard peak is a peak found in a system using a liquid chromatograph (LC) as the inlet, the title bar contains the following text:</p> <p>Peak Information – <i>component name</i></p> |

Table 24. Peak types for the Peak Information dialog box (Sheet 2 of 2)

| Peak type | Description |
|---|--|
| Qualifier ion peak (GC) | In a system using a gas chromatograph (GC) as the inlet, if the peak is for a Qualifier Ion, the title bar contains the following text: Peak Information – <i>component name</i> – Qual Ion Mass <i>xxx.x</i> , where <i>xxx.x</i> represents the mass of the qualifier ion |
| Spectrum candidate peak (GC) (First candidate) | In a system using a gas chromatograph (GC) as the inlet, if the peak is for the first spectrum candidate, the title bar contains the following text: Peak Information – <i>component name</i> – <i>spectrum candidate</i> , where <i>spectrum candidate</i> is the first candidate |
| Spectrum Candidate Peak (GC) (Second and Third Candidate) | In a system using a gas chromatograph (GC) as the inlet, if the peak is for the first spectrum candidate, the title bar contains the following text: Peak Information – <i>component name</i> – <i>spectrum candidate</i> , where the spectrum candidate is the second or third candidate |

Table 25. Peak Information dialog pages based on the peak type

| Peak Type | Info | Flags | More Flags | Suitability | Spectrum | More Info | Chro |
|---|-----------------|-------|------------|-------------|----------|-----------|------|
| No peak | – [*] | – | – | – | – | – | – |
| Standard peak (LC) | X ^{**} | X | X | X | X | – | – |
| Qualifier ion peak (GC) | X | X | X | X | X | X | X |
| Spectrum candidate peak (GC) (First candidate) | X | X | X | X | X | X | – |
| Spectrum Candidate Peak (GC) (Second and Third Candidate) | – | – | – | – | X | X | X |

^{*} (–) This page is not displayed.

^{**} (X) This page is displayed.

No Peak Page – Peak Information Dialog Box

If the Xcalibur data system does not find a peak for the component, only the No Peak page is displayed with the following message:

No Peak Found. Cannot show Peak Info.

Info Page – Peak Information Dialog Box

Use the Info page of the Peak Information dialog box to review the following peak properties for the current component.

This table describes the parameters on the Info page of the Peak Information dialog box.

Table 26. Info page parameters – Peak Information dialog box

| Parameter | Description |
|---------------------|---|
| Left (min) | Displays the left extreme of the integration baseline for the current component in minutes (read-only). |
| Apex (min) | Displays the apex point in minutes of the integration baseline for the current component (read-only). |
| Right (min) | Displays the right extreme of the integration baseline for the current component in minutes (read-only). |
| Height | Displays the height of the current component peak apex in units of counts (read-only). |
| Area (cts-sec) | Displays the area of the current component peak in units of count-seconds (read-only). |
| Baseline | Displays the baseline height directly below the apex of the current component peak in units of counts (read-only). |
| Base Peak (m/z) | Displays the mass-to-charge ratio of the ion with the largest response in the current component peak (read-only). |
| Signal To Noise | Displays the measured signal-to-noise ratio at the apex of the current component peak. (read-only) |
| Expected RT (min) | Displays the expected retention time of the current component (read-only). |
| ISTD Response | Displays the integrated area in units of count-minutes or the height of the apex in units of counts for the current component peak (read-only). |
| Response Ratio | Displays the ratio of the peak sample peak area or height to the internal standard area or height (read-only). |
| Calculated Amount | Displays the amount of sample calculated by the Xcalibur data system using the response ratio and the calibration curve (read-only). |

Flags Page – Peak Information Dialog Box

Use the Flags page of the Peak Information dialog box to review the following peak properties for the current component.

This table describes the parameters on the Flags page of the Peak Information dialog box.

Table 27. Flags page parameters – Peak Information dialog box (Sheet 1 of 4)

| Parameter | Description |
|-------------------------|--|
| Integration info | |
| Detected By | <p>Displays the method that the data system used to detect the peak:</p> <p>Spectrum: The application uses the user-defined mass/intensity pairs and applies a spectral-matching algorithm to find the peak that contains the closest match to the comparison spectrum. Up to 50 entries are allowed to define the comparison spectrum.</p> <p>Highest Peak: The application searches for the highest peak within the search window.</p> <p>Nearest RT: The application searches for the peak nearest to the expected retention time.</p> |
| Left Edge Type | <p>Displays how the data system detected the left baseline edge of the current peak. The data system displays one of the following peak baseline detection methods:</p> <p>Edge Type Reported by Xcalibur: Peak Criteria Met</p> <p>Baseline (B): The edge of the peak is at baseline level.</p> <p>Valley (V): The edge of the peak is in a peak valley.</p> <p>Manual (M): The edge of the peak has been adjusted manually.</p> <p>Stripe (S): The edge of the peak reached the Constrain Peak Height Percent specified in the method.</p> <p>Tail (T): The edge of the peak reached the Constrain Peak Height Trailing Factor limit before the Height Percent.</p> <p>Tilt (-): An error occurred before the data system could determine the edge of the peak.</p> <p>Unknown (?): An unknown error occurred.</p> |

Table 27. Flags page parameters – Peak Information dialog box (Sheet 2 of 4)

| Parameter | Description | | | | | | | | | | | | | | | |
|-----------------|--|------------|-----------|------------|----|----------|----------|----|----------|------|----|------|----------|----|--------|--------|
| | <p>View the edge type on many reports using a one-letter abbreviation for the left edge type and a one-letter abbreviation for the right edge type, as follows:</p> <table><tr><th>Peak ID</th><th>Left Edge</th><th>Right Edge</th></tr><tr><td>BB</td><td>Baseline</td><td>Baseline</td></tr><tr><td>BT</td><td>Baseline</td><td>Tail</td></tr><tr><td>-B</td><td>Tilt</td><td>Baseline</td></tr><tr><td>SS</td><td>Stripe</td><td>Stripe</td></tr></table> | Peak ID | Left Edge | Right Edge | BB | Baseline | Baseline | BT | Baseline | Tail | -B | Tilt | Baseline | SS | Stripe | Stripe |
| Peak ID | Left Edge | Right Edge | | | | | | | | | | | | | | |
| BB | Baseline | Baseline | | | | | | | | | | | | | | |
| BT | Baseline | Tail | | | | | | | | | | | | | | |
| -B | Tilt | Baseline | | | | | | | | | | | | | | |
| SS | Stripe | Stripe | | | | | | | | | | | | | | |
| Valid | Displays whether the data system successfully detected the peak indicated by selecting the Valid check box. | | | | | | | | | | | | | | | |
| Right Edge Type | <p>Displays how the data system detected the right baseline edge of the current peak. The application displays one of the following peak baseline detection methods:</p> <p>Edge Type Reported by Xcalibur</p> <p>Baseline (B): The edge of the peak is at baseline level.</p> <p>Valley (V): The edge of the peak is in a peak valley.</p> <p>Manual (M): The edge of the peak has been adjusted manually.</p> <p>Stripe (S): The edge of the peak reached the Constrain Peak Height Percent specified in the method.</p> <p>Tail (T): The edge of the peak reached the Constrain Peak Height Trailing Factor limit before the Height Percent.</p> <p>Tilt (-): An error occurred before the edge of the peak could be determined.</p> <p>Unknown (?): An unknown error occurred.</p> | | | | | | | | | | | | | | | |
| | <p>View the edge type on many reports using a one-letter abbreviation for the left edge type and a one-letter abbreviation for the right edge type, as follows:</p> <table><tr><th>Peak ID</th><th>Left Edge</th><th>Right Edge</th></tr><tr><td>BB</td><td>Baseline</td><td>Baseline</td></tr><tr><td>BT</td><td>Baseline</td><td>Tail</td></tr><tr><td>-B</td><td>Tilt</td><td>Baseline</td></tr><tr><td>SS</td><td>Stripe</td><td>Stripe</td></tr></table> | Peak ID | Left Edge | Right Edge | BB | Baseline | Baseline | BT | Baseline | Tail | -B | Tilt | Baseline | SS | Stripe | Stripe |
| Peak ID | Left Edge | Right Edge | | | | | | | | | | | | | | |
| BB | Baseline | Baseline | | | | | | | | | | | | | | |
| BT | Baseline | Tail | | | | | | | | | | | | | | |
| -B | Tilt | Baseline | | | | | | | | | | | | | | |
| SS | Stripe | Stripe | | | | | | | | | | | | | | |

Table 27. Flags page parameters – Peak Information dialog box (Sheet 3 of 4)


| Parameter | Description |
|-------------------|---|
| Flags | |
| Saturated | Displays whether any of the scans in the peak were saturated. The data system indicates one or more saturated scans were detected by displaying in the Saturated check box. |
| Calculated Amount | Displays the amount of sample, as calculated by the data system using the response ratio and the calibration curve. |
| Valley Detect | Displays whether valley detection was turned on in the processing method. The data system indicates that valley detection was activated if the Valley Detect check box is selected. |
| QC Failed | <p>Displays whether a sample failed a QC check.</p> <p>If the calculated amount is greater than the specified percentage difference from the expected amount, then a sample fails the QC test. For example, if the tolerance level is 10% and the expected amount is 100%, calculated amounts less than 90% or greater than 110% fail.</p> <p>The data system selects the check box to indicate that the sample type was QC and that it failed the QC test.</p> |
| RT Ref OK | <p>Displays whether the data system found the retention time reference component and whether it was used correctly by the processing method.</p> <p>If there was a retention time reference, then the check box indicates whether the retention time reference peak was found. The data system indicates it found the peak by selecting the check box. The application indicates it looked for the peak but did not find it by leaving the box blank.</p> <p>If there is no retention time reference, then the check box is selected because there is no correction to be made.</p> |
| Response OK | <p>Displays whether a response factor was calculated.</p> <p>The data system selects the check box to indicate that it found the peak and the peak's internal standard and correctly calculated the response ratio.</p> |

Table 27. Flags page parameters – Peak Information dialog box (Sheet 4 of 4)

| Parameter | Description |
|---------------|---|
| Response Low | Displays whether the calculated amount for the peak was less than the lowest specified standard amount of the component in the calibration curve. In this case the calculated amount has been determined by extrapolation from the lowest level. The data system selects the check box to indicate that the amount was calculated by extrapolation. If you force or include the origin, the application defines the lowest level to be 0.0. |
| Response High | Displays whether the calculated amount for the peak was greater than the highest specified standard amount of the component in the calibration curve. In this case, the calculated amount has been determined by extrapolation from the highest level. The data system selects the check box to indicate that the amount was calculated by extrapolation. |

More Flags Page – Peak Information Dialog Box

Use the More Flags page of the Peak Information dialog box to review the following properties for the current component.

A red check, , means that the parameter failed to meet the specified tolerance setting.

This table describes the parameters on the More Flags page of the Peak Information dialog box.

Table 28. More Flags page parameters – Peak Information dialog box (Sheet 1 of 2)

| Parameter | Description |
|---|---|
| Detection Thresholds | |
| When creating a processing method, set these thresholds in the Data Flags dialog box. | |
| Area | Set a specified area tolerance value to compare with the peak area and determine if the peak area is outside or within the tolerance. |
| Height | Set a specified height tolerance value to compare with the peak height and determine if the peak height is outside or within the tolerance. |

Table 28. More Flags page parameters – Peak Information dialog box (Sheet 2 of 2)

| Parameter | Description |
|---|--|
| Calibration and Quantitation Flag Thresholds | |
| When creating a processing method, set these thresholds in the Calibration and Quantitation Flags dialog box. You can view and change these settings in the Calibration Settings dialog box (see “ Calibration Settings Dialog Box ” on page 98). | |
| R-Squared | Set a specified coefficient tolerance value to compare with the coefficient of determination and determine if the coefficient of determination is outside or within the specified tolerance. |
| Detection Limit | Set a specified detection threshold tolerance value to compare with the concentration of the quantified peak and determine if the detection threshold is outside or within the specified tolerance. |
| Linearity Limit | Set a specified linearity threshold tolerance value to compare with the linearity threshold for the concentration of the quantified peak to determine if the linearity threshold is outside or within the specified tolerance. |
| Quantitation Limit | Set a specified quantitation threshold tolerance value to compare with the quantitation threshold for the concentration of the quantified peak and determine if the quantitation threshold is outside or within the specified tolerance. |
| Carry over Limit | Set a specified carry-over threshold tolerance value to compare with the carry-over threshold for the concentration of the quantified peak and determine if the carry-over threshold is outside or within the specified tolerance. |

Chro Page – Peak Information Dialog Box

The Chro page of the Peak Information dialog box is a read-only page that displays the single mass chromatogram of the current qualifier ion. The view centers on the retention time of the peak apex and has View Width specified in the Component Identification view of the Processing Setup window. The mass of the qualifier ion is displayed in the title bar Qual Ion Mass xxx.x. You cannot change this display.

This page is available for GC/MS data only.

Info or More Info Page – Peak Information Dialog Box

The Info or More Info page of the Peak Information dialog box is displayed if you have conducted a spectrum search.

- If the main component peak was detected using the spectrum search, then this page has the More Info tab and displays only the Spectrum box. In this case, the pages displayed are Info, More Info, Flags, More Flags, Suitability, and Spectrum.
- If peaks other than the main component were detected using the spectrum search (second or third candidates), then this page has the Info tab and displays both the Spectrum box and the Peak Info box. In this case, the pages displayed are Info, Chro, and Spectrum.

Use this page to review the following peak properties for the current search component.

This table describes the parameters on the Info and More Info pages of the Peak Information dialog box.

Table 29. Info or More Info page parameters – Peak Information dialog box (Sheet 1 of 2)

| Parameter | Description |
|------------------|---|
| Spectrum Results | Displays the calculated Forward, Reverse, and Match results for the found peak (read-only). This box appears on the More Info page for Candidate #1 of a spectrum search and on the Info page for Candidate #2 and Candidate #3 of a spectrum search. |
| Forward | Displays the forward threshold result for the current component (read-only). |
| Reverse | Displays the reverse threshold result for the current component (read-only). |
| Match | Displays the match threshold result for the current component (read-only). |
| Peak Info | Displays the calculated peak characteristics for Candidate #2 or Candidate #3 of a spectrum search (read-only). The Left, Apex, and Right peak boundaries in minutes, and the Area and Height of the peak are given in units of counts. This box on the More Info page does not appear for Candidate #1 of a spectrum search. This information is provided on the More page that appears for Candidate #1. |
| Left | Displays the left extreme of the integration baseline for the current component in minutes (read-only). |
| Apex | Displays the apex point in minutes of the integration baseline for the current component (read-only). |

Table 29. Info or More Info page parameters – Peak Information dialog box (Sheet 2 of 2)

| Parameter | Description |
|-----------|--|
| Right | Displays the right extreme of the integration baseline for the current component in minutes (read-only). |
| Area | Displays the area of the current component peak in units of count-seconds (read-only). |
| Height | Displays the height of the current component peak apex in units of counts (read-only). |

More Info Page – Peak Information Dialog Box

Use the More Info page of the Peak Information dialog box to review the results of the following peak tests for the current component.

This page is available for GC/MS data only.

This table describes the parameters on the More Info page of the Peak Information dialog box.

Table 30. More Info page parameters – Peak Information dialog box (Sheet 1 of 2)

| Parameter | Description |
|---------------------------|--|
| Ion Coelution Test | |
| Passed | Displays whether the qualifier ion displayed in the title bar “Qual Ion Mass xxx.x” passed or did not pass the Ion Coelution test. If the check box is not selected, no Ion Ratio test was performed and the Ion Ratio Test box is not displayed. If the check box is selected, the Ion Ratio test was performed and the Ion Ratio Test box on the Mass Info page is displayed at the bottom of the Mass Info page. |
| Ion Ratio Test | |
| Passed | Displays whether the qualifier ion displayed in the title bar “Qual Ion Mass xxx.x” passed or did not pass the Ion Ratio test. The Xcalibur data system does not display this box if the current qualifier ion did not pass the Ion Coelution test. |

Table 30. More Info page parameters – Peak Information dialog box (Sheet 2 of 2)

| Parameter | Description |
|------------------|---|
| Target Ratio% | Displays the calculated Target Ratio Percentage that Xcalibur calculated during the Ion Ratio test (read-only). The application does not display this box if the current qualifier ion did not pass the Ion Coelution test. |
| Absolute Window% | Displays the calculated Absolute Window Percentage that Xcalibur calculated during the Ion Ratio test (read-only). The application does not display this box if the current qualifier ion did not pass the Ion Coelution test. |

Spectrum Page – Peak Information Dialog Box

This read-only page displays the spectrum of the current peak at the apex retention time. You cannot make adjustments to this display.

Suitability Page – Peak Information Dialog Box

The Suitability page of the of the Peak Information dialog box displays the results of specific tests that might have been performed (as specified by the processing method) on the component peak to determine its suitability to be considered a valid peak. Each test is listed with the result of that test displayed to the right. There are three possible responses for each test: Passed, Failed, and Not Tested.

This table describes the parameters on the Suitability page of the Peak Information dialog box.

Table 31. Suitability page parameters – Peak Information dialog box (Sheet 1 of 2)

| Parameter | Description |
|--------------------------|--|
| Suitability Flags | |
| Symmetrical | Displays the results of the Symmetrical test as Passed, Failed, or Not Tested. This test indicates whether the peak is symmetrical about the apex. |
| Resolution | Displays the results of the Resolution test as Passed, Failed, or Not Tested. This test indicates whether multiple peaks are resolved. If neither peak baseline endpoint is valley detected, then the resolution passes. |
| Peak Width | Displays the results of the Peak Width test as Passed, Failed, or Not Tested. This test indicates whether the peak width is within specified limits. |
| Tailing | Displays the results of the Tailing test as Passed, Failed, or Not Tested. This test indicates whether the peak has tailing. |

Table 31. Suitability page parameters – Peak Information dialog box (Sheet 2 of 2)

| Parameter | Description |
|-----------------------|--|
| Column Overload | Displays the results of the Column Overload test as Passed, Failed, or Not Tested. This test indicates whether it is likely that the column was overloaded during acquisition. This test is based on an analysis of the baseline and peak shape. |
| Baseline Clipping | Displays the results of the Baseline Clipping test as Passed, Failed, or Not Tested. This test indicates whether the baseline is clipped outside the peak. This can occur if the chromatogram was started or terminated prematurely. |
| Signal-to-noise Ratio | Displays the results of the Signal-to-noise test as Passed, Failed, or Not Tested. This test indicates whether the minimum signal-to-noise ratio criteria are met. |
| Concave | Displays the results of the Concave test as Passed, Failed, or Not Tested. This test indicates whether the peak exhibits a concave depression (local minimum) due to noise. |
| Saturation | Displays the results of the Saturation test as Passed, Failed, or Not Tested. This test indicates whether the detector was saturated during acquisition. |

Quantitation Results Sorting Order Dialog Box

Use the Quantitation Results Sorting Order dialog box to set the sort order for the samples in the results grid view of the Quan Browser window. The sort order defines the priority for each parameter used in the sort.

For more information, see [“Changing the Sort Order”](#) on [page 32](#).

This table describes the parameters in the Quantitation Results Sorting Order dialog box.

Table 32. Quantitation Results Sorting Order dialog box parameters (Sheet 1 of 2)

| Parameter | Description |
|----------------|--|
| Sorting | |
| First Order | Base the first sort order of the results grid view on any of the following column headings or file properties: <none>, %Difference, %RSD, Area/Height, Area/Height Ratio, Exclude, File Name, Integration Type, Level Name, Peak Status, Sample ID, Sample Type, or Acquisition Date. By default, the first order sort is set to the acquisition date of the file. You can select and sort with any of these sort options, even if the corresponding column is not currently displayed. For example, you can sort by sample type, even if you have selected the Sample Name check box in the Result List Column Hiding dialog box. |

Table 32. Quantitation Results Sorting Order dialog box parameters (Sheet 2 of 2)

| Parameter | Description |
|--------------------------|--|
| Sort in descending order | Sort the first criterion in descending (reverse) order. If you do not select this check box, the sort is in ascending order. |
| Second Order | Base the second sort order of the results grid view on any of the following column headings or file properties: <none>, %Difference, %RSD, Area/Height, Area/Height Ratio, Exclude, File Name, Integration Type, Level Name, Peak Status, Sample ID, Sample Type, or Acquisition Date. You can select and sort with any of these sort options, even if the corresponding column is not currently displayed. For example, you can sort by sample type, even if you have selected the Sample Name check box in the Result List Column Hiding dialog box. |
| Sort in descending order | Sort the second criterion in descending (reverse) order. If you do not select this check box, the sort is in ascending order. |
| Third Order | Base the third sort order of the results grid view on any of the following column headings or file properties: <none>, %Difference, %RSD, Area/Height, Area/Height Ratio, Exclude, File Name, Integration Type, Level Name, Peak Status, Sample ID, Sample Type, or Acquisition Date. You can select and sort with any of these sort options, even if the corresponding column is not currently displayed. For example, you can sort by sample type, even if you have selected the Sample Name check box in the Result List Column Hiding dialog box. |
| Sort in descending order | Sort the third criterion in descending (reverse) order. If you do not select this check box, the sort is in ascending order. |
| Button | |
| Save As Default | Save your current selection of second and third sort orders as your default set. The data system uses these sort orders to display the results grid view until you use the Quantitation Results Setting Order dialog box to change your column sorting preferences. |

Reports Dialog Box

Use the Reports dialog box to generate reports on each sample row within the current bracket.

The Reports dialog box opens with a selection of Report templates and data files preloaded. These are obtained from the processing method that was previously defined and loaded along with the result file. This means that the values loaded might change as various brackets are selected.

Note You can use the standard report templates provided with the Xcalibur data system or you can create your own custom report templates. For information on how to create a report template, refer to the *XReport User Guide*.

For more information, see “[Generating Reports](#)” on [page 68](#).

This table describes the parameters in the Reports dialog box.

Table 33. Reports dialog box parameters (Sheet 1 of 5)

| Parameter | Description |
|-----------------------|---|
| Sample Reports | |
| Enabled | <p>Specifies whether the reports marked with Yes in their row are processed using the template that appears in the Report Template Name box. For example, if Yes appears in the QCs and Unknowns boxes and the Enable box in the row displays Yes, then the data system prints these sample reports. If the Enabled box in the row is clear, then the application does not print any sample reports.</p> <p>When you click an Enabled box, a check box control appears. If you select this check box and you click another cell, the application displays the word <i>Yes</i> to indicate that the report is enabled. If you do not select the check box, the cell remains blank when you click another cell.</p> |

Table 33. Reports dialog box parameters (Sheet 2 of 5)

| Parameter | Description |
|-----------|--|
| Stds | <p>Specifies whether the Standards (Stds) sample reports marked with Yes are processed using the report template that appears in the Report Template Name box of the same row.</p> <p>To print a Standard sample report, the data system must display Yes in the Enabled box in the same row. For example, if the Std box displays Yes and the Enable box in the same row displays Yes, then the application prints the Std sample report. However, if the Enabled box in the row is clear, then the application does not print the Std sample report.</p> <p>When you click a Stds box, a check box control appears. If you select this check box and you click another cell, the data system displays the word <i>Yes</i> to indicate that the report is enabled for Standard samples. If it is not selected, the cell remains blank when you click another cell.</p> |
| QCs | <p>Specifies whether the Quality Controls (QCs) sample reports marked with Yes are processed using the report template that appears in the Report Template Name box of the same row.</p> <p>To print a Quality Controls sample report, the data system must display Yes in the Enabled box in the same row. For example, if the Std box displays Yes and the Enable box in the same row displays Yes, then the data system prints the QCs sample report. However, if the Enabled box in the row is clear, then the application does not print the QCs sample report.</p> <p>When you select a QCs box, a check box control appears. If you select this check box and you click another cell, the data system displays the word <i>Yes</i> to indicate that the report is enabled for QC samples. If you do not select the check box, the cell remains blank when you click another cell.</p> |

Table 33. Reports dialog box parameters (Sheet 3 of 5)

| Parameter | Description |
|-----------|--|
| Unks | <p>Specifies whether the Unknowns (Unks) sample reports marked with Yes are processed using the report template that appears in the Report Template Name box of the same row.</p> <p>To print an Unknowns sample report, the data system must display Yes in the Enabled box in the same row. For example, if the Unks box displays Yes and the Enable box in the same row displays Yes, then the application prints the Unks sample report. However, if the Enabled box in the row is clear, then the application does not print the Unks sample report.</p> <p>When you select an Unks box, a check box control appears. If you select this check box and you click another cell, the data system displays the word <i>Yes</i> to indicate that the report is enabled for Unk samples. If you do not select the check box, the cell remains blank when you click another cell.</p> |
| Other | <p>Specifies whether the Other sample reports marked with Yes are processed using the report template that appears in the Report Template Name box of the same row.</p> <p>To print an Other sample report, the data system must display Yes in the Enabled box in the same row. For example, if the Other box displays Yes and the Enable box in the same row displays Yes, then the application prints the Other sample report. However, if the Enabled box in the row is clear, then the application does not print the Other sample report.</p> <p>When you select an Other box, a check box control appears. If you select this check box and you click another cell, the application displays the word <i>Yes</i> to indicate that the report is enabled for Other samples. If you do not select the check box, the cell remains blank when you click another cell.</p> |

Table 33. Reports dialog box parameters (Sheet 4 of 5)

| Parameter | Description |
|------------------------|--|
| Save As | <p>Specifies the file type for the sample report. These are the file type selections:</p> <ul style="list-style-type: none"> • None: print only, no exported file • Text: ASCII plain text file (TXT) • Doc: Microsoft Word document (DOC) • HTML: hypertext markup language file (HTML) • PDF: portable document format file (PDF) • RTF: rich text format file (RTF) • XLS: Excel spreadsheet (XLS) <p>The data system saves the exported file with the sample file name and the appropriate extension in the folder where the selected result files are stored.</p> |
| Report Template Name | <p>Specifies the report template to be used in processing the data and generating the reports indicated by Yes in each row.</p> <p>Double-clicking the grid cell opens the Open Report Template dialog box, where you select a template file. For information on how to create a report template, refer to the <i>XReport User Guide</i>.</p> |
| Summary Reports | |
| Enabled | <p>Specifies whether the summary reports are processed using the templates that appear in the Report Template Name boxes. For example, if the Enable box in the row displays Yes, then the data system prints the summary report defined by the report template in the row. If the Enabled box in the row is clear, then the application does not print this summary report.</p> <p>When you select an Enabled box, a check box control appears. If you select this check box and you click another cell, the application displays the word <i>Yes</i> to indicate that the report is enabled. If you do not select the check box, the cell remains blank when you click another cell.</p> |

Table 33. Reports dialog box parameters (Sheet 5 of 5)

| Parameter | Description |
|------------------------|---|
| Save As | <p>Specifies the file type for the summary report. These are the valid export file types:</p> <ul style="list-style-type: none"> • None: print only, no exported file • Text: ASCII plain text file (TXT) • Doc: Microsoft Word document (DOC) • HTML: hypertext markup language file (HTML) • PDF: portable document format file (PDF) • RTF: rich text format file (RTF) • XLS: Excel spreadsheet (XLS) <p>The data system saves the exported file with the sample file name and the appropriate extension in the folder where the selected result files are stored.</p> |
| Report Template Name | <p>Specifies the summary report template to be used in processing the summary report.</p> <p>Double-clicking the grid cell opens the Open Report Template dialog box, where you select a template file. For information on how to create a report template, refer to the <i>XReport User Guide</i>.</p> |
| Other Controls | |
| Include Sample Report | Select whether to print sample reports when you process the current data. This check box controls the printing of all of the sample reports defined in the Sample Reports box. |
| Include Summary Report | Select whether to print summary reports when you process the current data. This check box controls the printing of all of the summary reports defined in the Summary Reports box. |
| Select Samples | Opens the Select Report Samples dialog box, where you select the samples to be included in the reports. For more information, see “Select Report Samples Dialog Box” on page 128 . |
| Print Reports | Print the reports for the samples you have selected. |

Result List Column Hiding Dialog Box

The Result List Column Hiding dialog box contains the list of possible columns that can be viewed in the results grid. Selecting the check box to the left of the column name adds the column to the grid, and clearing the check box removes the column from the grid.

This table describes the parameters in the Result List Column Hiding dialog box.

Table 34. Result List Column dialog box parameters

| Parameter | Description |
|------------------|--|
| Selected Columns | <p>These settings indicate whether to display a column in the results grid. The following columns are available:</p> <ul style="list-style-type: none">• File Name• Sample Type• Sample Name• Integration Type• Area/Height• ISTD Area/Height• Area/Height Ratio• Specified Amount• Calculated Amount• Percent Difference• Percent RSD• Peak Status• Levels• Units• Retention Time• Sample ID• Exclude |

Select Level Dialog Box

Use the Select Level dialog box to select a level for a Standard or QC sample type in the results grid view of the Quan Browser window.

Note Whenever you change a QC, Blank, or Unknown sample type to a Standard sample type or change a Standard, Blank, or Unknown sample type to a QC sample type in the Quan Browser window results grid view, the Select Level dialog box opens so that you can select one of the available levels for the Standard or QC sample.

This table describes the parameters in the Select Level dialog box.

Table 35. Select Level dialog box parameters

| Parameter | Description |
|---------------------|--|
| Levels | <p>Select one of the available levels for the Standard or QC sample whenever you change a QC, Blank, or Unknown sample type to a Standard sample type or change a Standard, Blank, or Unknown sample type to a QC sample type.</p> <p>When you select a level from the Levels list, the read-only information in the lower list displays the other component names that are assigned to the selected level and the amount assigned to the selected level for each component.</p> |
| Levels Table | |
| Name | View the components that are assigned the same level as the one currently selected in the Levels list. |
| Amount | View the Component Amount assigned to the level currently selected in the Levels list for the component name selected in the Name column. |

Select Report Samples Dialog Box

The Select Report Samples dialog box displays the samples in the current bracket or group in the Sample Choices box.

From the Sample Choices dialog box, you can pick which samples are to be processed when sample reports are selected. Use the SHIFT and CTRL keys to select multiple samples. The data system remembers the selected samples for each bracket until the application is terminated or a new file is opened.

This table describes the parameters in the Select Report Samples dialog box.

Table 36. Select Report Samples dialog box parameters

| Parameter | Description |
|-------------------------|--|
| Sample Choices | |
| Raw File | View or change the raw files in the current bracket or current group of samples. These are the sample files that you select from for processing and for printing a report. |
| Sample Type | View the sample type of the raw file displayed to the left on the same row. |
| Selected Samples | |
| Raw File | View or change the raw files that have been selected from the Sample Choices box. These are the sample files to be processed so that you can print a report. |
| Sample Type | View the sample type of the raw file displayed to the left on the same row. |
| Buttons | |
| Add | Add files selected in the Sample Choices box to the Selected Samples box. |
| Remove | Return files selected in the Selected Choices box back to the Sample Choices box. |
| Add All | Add all of the files in the Sample Choices box to the Selected Samples box. |
| Remove All | Return all files in the Selected Choices box back to the Sample Choices box. |

User Identification Settings Dialog Box

Use the User Identification Settings dialog box to select and test mass, scan filter, relative peak height threshold, peak identification, and peak integration settings. If these standard options do not provide the desired results, Quan Browser also provides advanced options using the Advanced page. The pages that are displayed depend on whether you are currently using the Genesis, the ICIS, or the Avalon peak detection algorithm.

For information about the identification, detection, integration, advanced, and flag parameters, see these topics:

- “Identification Page – User Identification Settings Dialog Box” on page 130
- “Detection Page – User Identification Settings” on page 134
- “Integration Page – User Identification Settings Dialog Box” on page 139
 - “Genesis Integration Page Parameters” on page 139
 - “ICIS Integration Page Parameters” on page 141
 - “Avalon Integration Page Parameters” on page 142
- “Advanced Page – User Identification Settings Dialog Box” on page 145
 - “Genesis Advanced Page Parameters” on page 145
 - “ICIS Advanced Page Parameters” on page 147
- “Flags Page – User Identification Settings Dialog Box” on page 148

❖ To open the User Identification Settings dialog box

Right-click the chromatogram view in the Quan Browser window and choose **User Peak Detection Settings** from the shortcut menu.

Identification Page – User Identification Settings Dialog Box

Use the Identification page to change the current component name, mass range, scan filter, and retention time for the selected component. You can then test the results of the new settings by clicking Apply or OK. You can apply the new settings to all files in the results grid by clicking Apply To All.

The Identification page helps to narrow the search parameters and to set filters so that the peak detection algorithms have an easier time of locating the peaks. The compound of interest is displayed in the Name box. This is a read-only field. To select a different compound, select the name in the component list on the right side of the display.

This table describes the parameters on the Identification page of the User Identification Settings dialog box.

Table 37. Identification page parameters – User Identification Settings dialog box (Sheet 1 of 5)

| Parameter | Description | | |
|-----------|--|----------------|------------------------|
| Name | Displays the component name selected in the components pane for the active processing method. | | |
| Plot Type | These three lists display the type of trace and optional trace math operation that is stored in the processing method. Only certain combinations of plot types are possible as shown in the following table: | | |
| | 1st Plot Type | Math Operation | 2nd Plot Type |
| | Mass Range | None | n/a |
| | Mass Range | ± | Mass Range |
| | TIC | | n/a |
| | TIC- | | Mass Range |
| | TIC- | | Base Peak |
| | Base Peak | None | n/a |
| | Base Peak | ± | Mass Range |
| | Analog 1, 2, 3, or 4 | None | n/a |
| | Analog 1, 2, 3, or 4 | ± | Digital* 1, 2, 3, or 4 |
| | Digital 1, 2, 3, or 4 | None | n/a |
| | Digital 1, 2, 3, or 4 | ± | Digital* 1, 2, 3, or 4 |

Table 37. Identification page parameters – User Identification Settings dialog box (Sheet 2 of 5)

| Parameter | Description |
|-------------|---|
| Scan Filter | <p>This box displays the current scan filter for the active raw file (RAW). You can use a scan filter to specify that processing is to be applied to a subset of the scans in a raw file.</p> <p>To apply a different scan filter, select a new filter from the scan filter list (most common method), select a new filter from the list and edit the scan filter, or type a new scan filter command string into the box using the scan filter format.</p> <p>For example, the following scan filter:</p> <p style="padding-left: 40px;">c full ms [26.81–251]</p> <p>finds all scans in a raw file that have the following properties:</p> <p style="padding-left: 40px;">centroid data</p> <p style="padding-left: 40px;">Scan Mode: Full</p> <p style="padding-left: 40px;">Scan Power: MS</p> <p style="padding-left: 40px;">Product Ion Mass Range: <i>m/z</i> 26.81 to 251.00</p> <p>The data system displays this box when you select a Base Peak trace for an MS detector type. The box displays the range within which the application is to search for the highest peak.</p> <p>If you type a single <i>m/z</i> value in this box, that defines the base peak.</p> <p>To change the base peak mass range, type the value in the box. A mass range from <i>m/z</i>=A to <i>m/z</i>=B is entered in the format A–B.</p> <p>The data system displays this box when you select a Base Peak ± Mass Range trace combination for an MS detector type. The box displays the mass range for the second, Mass Range, trace type.</p> <p>To change the range or to add a new range, type the range in the box. The format is <i>Low Mass–High Mass</i>. For example, for the range <i>m/z</i> 123 through 456, type: 123–456.</p> |

Table 37. Identification page parameters – User Identification Settings dialog box (Sheet 3 of 5)

| Parameter | Description |
|---|---|
| Mass | Displays the masses stored in the processing method. This display area changes to accommodate the type of data required. When a single mass range is required, there is a single edit box displaying the current value. If two mass ranges are required (such as the case of a trace defined as a Mass Range \pm Mass Range or Base Peak \pm Mass Range), this box is replaced by two boxes (in the case of Base Peak \pm Mass Range, this box is replaced by the BP and MR boxes). In the case of a TIC (no trace operator in use), Analog, or Digital traces, this box is blank. |
| Keys (read-only) | Displays user comments about the processing method. |
| Retention Time | |
| The settings in this area define the expected retention time in minutes of the component peak and the error window in seconds for the retention time. | |
| Expected | <p>This box displays the expected peak width parameter (in seconds). This controls the minimum width that a peak is expected to have if valley detection is enabled.</p> <p>With valley detection enabled, any valley points nearer than the <i>expected width/2</i> to the top of the peak are ignored. If a valley point is found outside the expected peak width, the data system terminates the peak at that point. It always terminates a peak when the signal reaches the baseline, independent of the value set for the expected peak width. The valid range is 0.0 to 999.0 seconds. To change the current value, type a new width in the Expected box.</p> |
| Window | This box displays the allowable retention time window for the elution of the selected component. The valid range is 1.0 to 999.0 seconds. To change the time window or to enter a new time window, type the number of seconds in the (retention time) Window box. |
| Use as RT Reference (read-only) | This check box indicates whether the actual retention time (RT) of the active component [as displayed in the Name combo box in the same view] was used to adjust the expected retention time of another component. |
| View Width | This box displays the current view width (in minutes). The valid range is dependent upon the configured hardware. To change the view width, type the desired time in the View Width text box. |

Table 37. Identification page parameters – User Identification Settings dialog box (Sheet 4 of 5)

| Parameter | Description |
|-----------------------------------|--|
| Adjust Using (read-only) | This check box indicates whether the expected retention time (RT) of the active component (as displayed in the Name box in the same view) is to be adjusted using the actual retention time of the RT Reference, such as an internal standard. The data system displays the RT Reference in the Adjust Using box to the right of this check box. |
| Adjust Using (box) (read-only) | This box displays the retention reference component that data system uses to adjust the expected retention time of the active component (as displayed in the Name box in the same view. The data system uses the actual retention time of the RT Reference component to correct the retention time of the active component. The application provides the following correction to the expected retention time. $\text{Comp RT}_{\text{adjusted}} = \frac{\text{Comp RT}_{\text{expected}} \times \text{Ref RT}_{\text{actual}}}{\text{Ref RT}_{\text{expected}}}$ <p>Where:</p> <p>Comp RT_{adjusted} is the adjusted expected RT of the active component.</p> <p>Ref RT_{actual} is the actual RT of the reference component.</p> <p>Ref RT_{expected} is the expected RT of the reference component.</p> |
| Detector Type | |
| Type | View detector type options. These are the detectors that you have configured using the Foundation Instrument Configuration dialog box. |
| Peak Detection Algorithm | This list contains three options from which you can select a peak detection algorithm to recalculate the data using that algorithm. |

Table 37. Identification page parameters – User Identification Settings dialog box (Sheet 5 of 5)

| Parameter | Description |
|----------------|---|
| Buttons | |
| Apply | Applies the current peak detection parameters to the selected component of the selected sample in the current sequence. |
| Apply to All | Applies the current peak detection parameters to all samples that are currently displayed in the sequence. For example, if “Standards” are displayed, the current peak detection parameters are applied to only the Standards samples and not to any other samples. If “All” samples are displayed, the current peak detection parameters are applied to all samples in the current sequence. If one or more Standard samples are changed, the data system recalculates all quantitation parameters, including peak areas and the calibration curve. If one or more Standard samples are changed, the data system recalculates all quantitation parameters, including peak areas and the calibration curve. |

Detection Page – User Identification Settings

Use the Detection page from the User Identification Settings dialog box to change the current peak detection criteria for component detection. The parameters on the Detection page vary depending on whether you use a GC or LC, whether the detection method is Spectrum, Highest Peak, or Nearest RT, and the selected integration algorithm.

This table describes the parameters on the User Identification Settings – Detection page.

Table 38. Detection page parameters – User Identification Settings dialog box (Sheet 1 of 5)

| Parameter | Description |
|---------------------|---|
| Highest Peak | This option is available for both LC/MS and GC/MS data. Use the highest peak in the chromatogram for component identification. |
| Nearest RT | This option is available for both LC/MS and GC/MS data. Use the peak with the nearest retention time in the chromatogram for component identification. |
| Minimum Peak Height | This parameter is available for the ICIS and Genesis algorithms. View or change the peak signal-to-noise criteria that needs to be equaled or exceeded for the data system to use the Nearest RT Peak Identification criteria. When identifying components, the application ignores all chromatogram peaks that have signal-to-noise values that are less than the S/N Threshold value. The valid range is 0.0 (all peaks) through 999.0. |

Table 38. Detection page parameters – User Identification Settings dialog box (Sheet 2 of 5)

| Parameter | Description |
|---|--|
| Spectrum | <p>This option is only available for GC/MS data. You can use a reference spectrum defined in the processing method for component identification. The data system attempts to match the reference spectrum with a series of unknown spectra and calculates a score value for each comparison.</p> <p>If you select the Spectrum option for GC/MS data, the following parameters appear:</p> <ul style="list-style-type: none"> • Spectrum Peak Detection: Displayed only for GC/MS data when you select the Spectrum option. You must also have selected the MS detector type on the Identification page. • Spectrum Peak Identification table: Enter mass/charge [m/z] and intensity data for up to 50 spectrum peaks. The data system uses this data to identify the active component in the Find algorithm. It displays this table only when you select the Spectrum option for GC/MS data. • m/z: View the mass/charge [m/z] value for one peak in the reference spectrum. The intensity for this m/z value is given in the adjacent Intensity table box. Use the other rows of the table to enter as many as 50 m/z values. The data system uses this data to identify the active component in the Find algorithm. It displays this table only when you select the Spectrum option for GC/MS data. • Intensity: Enter intensity data for one peak in the reference spectrum. The m/z value for this intensity is given in the adjacent m/z Table box. Use the other rows of the table to enter as many as 50 intensity values. The data system uses this data to identify the active component in the Find algorithm. It displays this table only when you select the Spectrum option for GC/MS data. |
| Thresholds (for the Spectrum option) | |
| Forward | Set a threshold value for Forward comparisons between the reference spectrum and candidates in the chromatogram. A Forward search is a direct matching algorithm comparing unknowns against the reference spectrum in the peak identification table. The match is scored on a scale of 0 to 999. A perfect match results in a score of 999. As a general guide, 900 or greater is an excellent match; 800 to 900 a good match; 700 to 800 a fair match. Less than 600 is a poor match. Unknown spectra with many peaks tend to score lower than similar spectra with fewer peaks. |
| Reverse | Set a threshold value for Reverse comparisons between the reference spectrum and candidates in the chromatogram. A Reverse search ignores any peaks in the unknown that are not in the reference spectrum in the peak identification table. The match is scored on a scale of 0 to 999. A perfect match results in a score of 999. As a general guide, 900 or greater is an excellent match; 800 to 900 a good match; 700 to 800 a fair match. Less than 600 is a poor match. A spectrum with many peaks tends to score more highly in a Reverse match than a Forward match. |

Table 38. Detection page parameters – User Identification Settings dialog box (Sheet 3 of 5)

| Parameter | Description |
|---|---|
| Match | Set a threshold value for Match comparisons between the reference spectrum and candidates in the chromatogram. Match is scored on a scale of 0 to 999. The Match algorithm is a complex probability factor based on the differences between the Forward factors of all the candidates. If one candidate has a Forward matching factor of 900 and the next best is only 300, the probability of the component being correctly identified is high and so the Match factor is scored highly for the first candidate. If the Forward factors for all the candidates are similar, whether high or low, the Match factor is low. |
| Ion Ratio Confirmation (GC/MS data) | |
| If you are working with GC/MS data and you select the Highest Peak or Nearest RT options, the Ion Ratio Confirmation area appears. You must also have selected the MS detector type on the Identification page. | |
| Enable | Indicates whether Ion Ratio Confirmation is enabled. |
| Ion Ratio Using: Area or Height | View the currently selected peak quantitation method: area or height. The data system uses the same method to calculate qualifier ion peak response and then target ratio. You can change this parameter by selecting the Area or Height options in the Response box on the Calibration page. |
| Qualifier Ion Table | <p>Specify up to five qualifier ions in this box to confirm the detection of a target analyte. You can also set the coelution window and select a method for calculating the target ion ratio window and tolerance.</p> <p>Use this table to enter mass/charge [m/z] and target ratio tolerances [Window ($\pm\%$)] data for up to five qualifier ions.</p> <ul style="list-style-type: none"> • If you are using Area response, the data system integrates each qualifier ion peak and ratios it with the quantitation peak area. The application then compares this ratio with your specified target ratio. If the calculated ratio is outside of the target ratio by more than your specified tolerance [Window ($\pm\%$)], the quantitation peak is rejected. • If you are using Height response, the data system ratios the qualifier ion peak height with that of quantitation peak. The application then compares this ratio with your specified target ratio. If the calculated ratio is outside of the target ratio by more than your specified tolerance [Window ($\pm\%$)], the quantitation peak is rejected. |
| m/z column | This column in the Qualifier Ion table displays the mass/charge [m/z] value for a qualifier ion. The target ratio tolerance for this m/z value is displayed in the adjacent Window($\pm\%$) table box. Use other rows of the table to enter data for five qualifier ions. The data system uses this data to confirm the identity of a quantitation peak by comparing the relative responses of qualifier ion and quantitation peaks with predetermined values. |

Table 38. Detection page parameters – User Identification Settings dialog box (Sheet 4 of 5)

| Parameter | Description |
|--------------|---|
| Target Ratio | This column in the Qualifier Ion table displays the Target Ratio (%) value for a qualifier ion. The m/z value and target ratio tolerance for the qualifier ion are given in the adjacent m/z and [Window ($\pm\%$)] table boxes. Use other rows of the table to enter data for five qualifier ions. The data system uses this data to confirm the identity of a quantitation peak by comparing the relative responses of qualifier ion and quantitation peaks with predetermined values. |
| Window | Use this column in the Qualifier Ion table to specify the Target Ratio tolerance for a qualifier ion. Use other rows of the table to enter data for five qualifier ions. The data system uses this data to confirm the identity of a quantitation peak by comparing the relative responses of qualifier ion and quantitation peaks with the specified values. |
| Window (%) | <p>Select Relative or Absolute.</p> <ul style="list-style-type: none"> Relative <p>Specify that the target ratio tolerance values in the Window ($\pm\%$) column of the qualifier ion table are relative values.</p> <p>For example, if the target ratio is 50% and the Window ($\pm\%$) parameter is 20%, the expected target ion ratio range is 40% to 60% (with the Absolute option this would be 30% to 70%). If the ion ratio is outside this range, the ion ratio confirmation test has failed and the data system sets the IRC Flag to false. If the qualifier ion peak/quantitation peak ratio is within range, the ion ratio confirmation test passes and the application sets the IRC Flag to true. The response of all specified qualifier ions must be within the respective ratio ranges for IRC to succeed.</p> <p>In assessing a target ion ratio range, the application truncates the range at 0% to avoid negative values.</p> Absolute <p>Specify that the target ratio tolerance values in the Window ($\pm\%$) column of the qualifier ion table are absolute values.</p> <p>For example, if the target ratio is 50% and the Window ($\pm\%$) parameter is 20%, the expected target ion ratio range is 30% to 70% (with the Relative option this would be 40% to 60%). If the qualifier ion peak/quantitation peak ratio is outside this range, the ion ratio confirmation test has failed and the data system sets the IRC Flag to false. If the qualifier ion peak/quantitation peak ratio is within range, the ion ratio confirmation test passes and the application sets the IRC Flag to true. The response of all specified qualifier ions must be within the respective ratio ranges for IRC to succeed.</p> <p>In assessing a target ion ratio range, the data system truncates the range at 0% to avoid negative values.</p> |

Table 38. Detection page parameters – User Identification Settings dialog box (Sheet 5 of 5)

| Parameter | Description |
|---------------------|---|
| Qualifier Coelution | <p>Qualifier Ion Coelution window.</p> <ul style="list-style-type: none"> • Prior to ion ratio confirmation, the data system generates a mass chromatogram for each specified qualifier ion. Each of these chromatograms must feature a peak, matching that of the quantitation masses. If the retention time of the qualifier ion peak apex lies outside of the Qualifier Ion Coelution window (centered on the quantitation peak), the application rejects the quantitation peak. • Quantitation peaks with matching qualifier ion peaks (within the Coelution window) are tested by Xcalibur for ion ratio confirmation. |
| Buttons | |
| Apply | <p>Apply the current peak detection parameters to the selected component of the selected sample in the current sequence.</p> <p>If one or more standard samples are changed, the data system recalculates all quantitation parameters, including peak areas and the calibration curve.</p> |
| Apply To All | <p>Apply the current peak detection parameters to all samples that are currently displayed in the sequence. For example, if you select the Standards option, the data system applies the current peak detection parameters only to the standards samples and not to any other samples. If you select the All samples option, the application applies the current peak detection parameters to all samples in the current sequence.</p> <p>If one or more standard samples are changed, the data system recalculates all quantitation parameters, including peak areas and the calibration curve.</p> |

Integration Page – User Identification Settings Dialog Box

Use the Integration page to change the current peak integration criteria for the selected component. You can then test the results of the new criteria by clicking Apply or OK. You can apply the new criteria to all files in the Result list by clicking Apply To All. These settings are used by the Detection algorithm.

Depending on the peak detection algorithm that you are using, one of three Integration pages is available. For information about the parameters on these pages, see these topics:

- [Genesis Integration Page Parameters](#)
- [ICIS Integration Page Parameters](#)
- [Avalon Integration Page Parameters](#)

Genesis Integration Page Parameters

This table describes the parameters on the Genesis Integration page of the User Identification Settings dialog box.

Table 39. Genesis Integration page parameters (Sheet 1 of 2)

| Parameter | Description |
|--------------------------|--|
| Smoothing Points | Determine the degree of data smoothing to be performed on the active component peak prior to peak detection and integration. The valid range is any odd value between 1 (no smoothing) through 15 (maximum smoothing). To smooth your component peak data prior to integration, type a value in the Smoothing Points box. See also Avalon Integration Page Parameters and ICIS Integration Page Parameters . |
| S/N Threshold | View or change the current signal-to-noise threshold for peak integration. Peaks with signal-to-noise less than this value are not integrated. Peaks with signal-to-noise greater than this value are integrated. The valid range is 0.0 to 999.0. To change the current value, type a new value in the S/N Threshold box. |
| Valley Detection Enabled | Use the Xcalibur 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. To turn this method on, select the Valley Detection check box. To turn this method off, ensure that the check box is clear. |

Table 39. Genesis Integration page parameters (Sheet 2 of 2)

| Parameter | Description |
|----------------------|---|
| Expected Width (sec) | <p>View the expected peak width parameter (in seconds). This controls the minimum width that a peak is expected to have if valley detection is enabled.</p> <p>With valley detection enabled, any valley points nearer than the <i>expected width/2</i> to the top of the peak are ignored. If a valley point is found outside the expected peak width, the data system 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. The valid range is 0.0 to 999.0 seconds. To change the current value, type a new width in the Expected Width box.</p> |
| Constrain Peak Width | <p>Limit 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. To constrain a peak width, select the Constrain Peak Width check box. The Peak Ht (%) box and the Tailing Factor box become available.</p> |
| Peak Ht (%) | <p>View or adjust the percent of the total peak height (100%) that a signal needs to be above the baseline before integration is turned on or off. This box is active only when the Constrain Peak Width check box is selected. The valid range is 0.0 to 100.0%. To enter this height, type the appropriate value in the Peak Ht (%) box.</p> |
| Tailing Factor | <p>View or adjust a factor that controls how the Xcalibur data system integrates the tail of a peak. This tailing factor is the maximum ratio of the trailing edge to the leading side of a constrained peak. This box is active only when the Constrain Peak Width box is selected. The valid range is 0.5 through 9.0.</p> |

ICIS Integration Page Parameters

This table describes the ICIS Integration page parameters.

Table 40. ICIS Integration page parameters – User Identification Settings dialog box

| Parameter | Description |
|----------------------|--|
| Smoothing Points | Type the number of points used in the moving average used to smooth the data. The valid range is any odd value from 1 through 15 points. The default value is 1 point. This value is used by the ICIS peak detection algorithm. |
| Baseline Window | Specify the number of scans over which to look for a local minima. The valid range is 1 through 500. The default value is 40 scans. This value is used by the ICIS peak detection algorithm. |
| Area Noise Factor | Specify the noise level multiplier used to determine the peak edge after the location of the possible peak. The valid multiplier range is 1 through 500. The default multiplier is 5. This value is used by the ICIS peak detection algorithm. |
| Peak Noise Factor | Specify the noise level multiplier used to determine the potential peak signal threshold. The valid multiplier range is 1 through 1000. The default multiplier is 10. This value is used by the ICIS peak detection algorithm. |
| Constrain Peak Width | Limit 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. To constrain a peak width, select the Constrain Peak Width check box. The Peak Ht (%) box and the Tailing Factor box become available. |
| Peak Ht (%) | View or adjust the percent of the total peak height (100%) that a signal needs to be above the baseline before integration is turned on or off. This box is active only when the Constrain Peak Width check box is selected. The valid range is 0.0 to 100.0%. To enter this height, type the appropriate value in the Peak Ht (%) box. |
| Tailing Factor | View or adjust a factor that controls how the Xcalibur data system integrates the tail of a peak. This tailing factor is the maximum ratio of the trailing edge to the leading side of a constrained peak. This box is active only when the Constrain Peak Width check box is selected. The valid range is 0.5 through 9.0. |

Avalon Integration Page Parameters

This table describes the parameters on the Avalon Integration page of the User Identification Settings dialog box.

Table 41. Avalon Integration page parameters (Sheet 1 of 3)

| Parameter | Description |
|--------------------------|--|
| Auto Calc Initial Events | <p>This button is active with the event list of the Avalon peak detection algorithm only if a raw file is open. When you click the button, Avalon automatically estimates the initial values for the detection of peaks based on the data in the current raw file, and then displays those initial values in the event list. Use this button to force Avalon to search for the best values of initial events that detect peaks in the data. Any timed event in the event list is unchanged when you click this button.</p> <p>Auto Calculate Initial Events determines initial values for the following events only: Start Threshold, End Threshold, Area Threshold, P-P [Resolution] Threshold, Bunch Factor, Negative Peaks, and Tension. Additionally, the user can specify timed events for these events in the same event list.</p> |
| Smoothing Points | View or adjust the number of points that the Xcalibur data system uses for chromatogram smoothing. The valid range for smoothing points is from 3 to 15. The number of smoothing points must be odd. To change the number of smoothing points, type the new number of points in the Smoothing Points box. |
| Event List | |
| Event List | <p>To detect peaks, Avalon uses the settings for initial events and user-defined timed events in the event list. To calculate values for initial events, click Auto Calc Initial Events.</p> <p>The event list in the Avalon Event List dialog box contains two hidden columns of information that are used by the algorithm and cannot be changed by the user: Event OP Code and Value2.</p> <p>There are seven initial entry integration events, which are identified by the initial value setting in the Time column. These are the default integration events required by the Avalon integration algorithm. You can change the Value of an initial entry integration event, but you cannot delete it or change its time value.</p> |
| Time | This column contains either the term initial value or a value of time in minutes. |

Table 41. Avalon Integration page parameters (Sheet 2 of 3)

| Parameter | Description |
|-------------------------|---|
| Event | View descriptions of detection parameters for initial events and timed events. You cannot change an event associated with an initial value. |
| Value | View the values associated with initial events or timed events. The range of factors allowed for each value is specific to each event. |
| Event List entry | |
| Time | View or change the currently highlighted entry from the Time column in the event list. |
| Event | View the currently highlighted entry in the Event column of the event list. An event cannot be changed that is listed with an Initial Value in the Time column. The Threshold and Bunch Factor parameters are the most important ones in controlling peak detection. |
| Start/End Threshold | Directly related to the RMS noise in the chromatogram, this is Threshold, the fundamental control used for peak detection. |
| Bunch Factor | The Bunch Factor is the number of points grouped together during peak detection. It controls the bunching of chromatographic points during integration and does not affect the final area calculation of the peak. The Bunch Factor must be an integer between 1 and 6; a high bunch factor groups peaks into clusters. |
| Area Threshold | Controls the area cutoff. Any peaks with a final area less than the area threshold is not detected. This control is in units of area for the data. |
| 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. This is specified as a percent of peak height overlap. |
| Negative Peaks | Automatically resets after a negative peak has been found. |
| Tension | Controls how closely the baseline should follow the overall shape of the chromatogram. A lower tension traces the baseline to follow changes in the chromatogram more closely. A high baseline tension follows the baseline less closely, over longer time intervals. Set in minutes. |

Table 41. Avalon Integration page parameters (Sheet 3 of 3)

| Parameter | Description |
|---------------------|---|
| 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 (solvent). 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. |
| Shoulders On | Turns on the detection of shoulders. |
| Shoulders Off | Shoulders Off Turns off the detection of shoulders. |
| Force Cluster On | Turns on the grouping of peaks into a single peak. |
| Force Cluster Off | Turns off the grouping of peaks into a single peak. |
| Disable Cluster On | Enables the grouping effect in the specified time range. |
| Disable Cluster Off | Disables the grouping effect in the specified time range. |
| Value | View or change the currently highlighted entry from the Value column in the event list. The range of factors allowed for each value is specific to each event. |
| Page Buttons | |
| Add | Add to add a time/event/value entry for a timed event in the event list. When you click Add, both the event list and the chromatogram display update automatically with the added specification in the currently selected chromatogram. |
| Delete | Delete to remove a highlighted event from the event list. You cannot delete initial values. |
| Change | Change to update a highlighted time/event/value entry in the event list. When you click Change, both the event list and the chromatogram display update automatically with the revised specification in the currently selected chromatogram. For initial events, only the values (and not the events) can be changed. |

Advanced Page – User Identification Settings Dialog Box

Use the Advanced page of the User Identification Settings dialog box to change the current advanced component detection criteria. Use these additional criteria if the standard detection criteria do not provide the desired results. You can then test the results of the new criteria by clicking Apply or OK. You can apply the new criteria to all files in the Result list by clicking Apply To All.

Advanced parameters used for the detection and integration of peaks are less often used but can provide adequate peak detection with the default parameters.

You can set advanced parameters for the Genesis and ICIS algorithms. For more information, see these topics:

- [Genesis Advanced Page Parameters](#)
- [ICIS Advanced Page Parameters](#)

Genesis Advanced Page Parameters

This table describes the parameters on the Genesis Advanced page of the User Identification Settings dialog box.

Table 42. Genesis Advanced page parameters (Sheet 1 of 2)



| Parameter | Description |
|------------------------|--|
| Report Noise As | |
| RMS | Select this option to calculate noise as RMS. |
| Peak To Peak | Select this option to calculate noise as peak-to-peak. |
| Manual Noise Region | Specify the region of the chromatogram that the Xcalibur data system uses to determine noise. You can click the Manual Noise Region icon,  , and drag the cursor horizontally across the region of the chromatogram that you want to select as the noise region or type the retention time (RT) in the RT Range box. The data system marks the region with a red baseline. |
| RT Range | Specify the retention time (RT) range. The RT range should be within the chromatogram range. You can click the Manual Noise Region icon,  , and drag the cursor horizontally across the region of the chromatogram that you want to select as the noise region or type a value in the RT Range box. The data system marks the region with a red baseline. |

Table 42. Genesis Advanced page parameters (Sheet 2 of 2)

| Parameter | Description |
|---------------------------------|--|
| Rise Percentage | View or adjust the percentage that the peak trace can rise above the baseline after passing through a minimum (before or after the peak). If the trace exceeds this value, the data system applies valley detection peak integration criteria. This test is applied to both the left and right edge of the peak. This criteria is useful for integrating peaks with long tails. The valid range is 0.1 to 500.0. To change the rise percentage, type a value in the Rise Percentage box. Click OK to apply the new peak detection criteria. |
| Valley S/N | View or adjust the signal-to-noise criteria that the data system uses for valley detection. The valid range is 1.0 to 100.0. To change the valley detection signal-to-noise criteria, type a value in the Valley S/N box. Click OK to apply the new peak detection criteria. |
| Peaks S/N Cutoff | View or adjust the signal-to-noise below which the data system defines the peak edge. For example, if the signal-to-noise at the apex is 500 and the Peak S/N Cutoff value is 200, the application defines the right and left edges of the peak when the S/N reaches a value less than 200. The valid range is 50.0 to 10 000.0. |
| Baseline Noise Tolerance (%) | View or adjust a value that controls how the baseline is drawn in the noise data. The higher the baseline noise tolerance value, the higher the baseline is drawn through the noise data. The valid range is 0.0 to 100.0. |
| Min Number Of Scans In Baseline | View or adjust the minimum number of scans that the Xcalibur data system uses to calculate a baseline. A larger number includes more data in determining an averaged baseline. The valid range is 2 to 100.0. |
| Number Of Background Scans | View or adjust the number of background scans used to determine the background. The valid range is 1 to 100. |

ICIS Advanced Page Parameters

This table describes the parameters on the ICIS Advanced page of the User Identification Settings dialog box.

Table 43. ICIS Advanced page parameters (Sheet 1 of 2)



| Parameter | Description |
|----------------------|---|
| Manual Noise Region | <p>Specify the region of the chromatogram that the Xcalibur data system uses to determine noise.</p> <p>When you select this check box, the RT Range box becomes available.</p> <p>You can click the Manual Noise Region icon, , and drag the cursor horizontally across the region of the chromatogram that you want to select as the noise region or type the retention time (RT) in the RT Range box. The data system marks the region with a red baseline.</p> |
| RT Range | <p>Specify the retention time (RT) range. The RT range should be within the chromatogram range.</p> <p>You can click the Manual Noise Region icon, , and drag the cursor horizontally across the region of the chromatogram that you want to select as the noise region or type a value in the RT Range box. The data system marks the region with a red baseline.</p> |
| Noise Method | |
| INCOS Noise | Use a single pass algorithm to determine the noise level. This value is used by the ICIS peak detection algorithm. |
| Repetitive Noise | Use a multiple pass algorithm to determine the noise level. This value is used by the ICIS peak detection algorithm. In general, this algorithm is more accurate in analyzing the noise than the INCOS noise algorithm, but it takes longer. |
| RMS | Specify that the data system calculates noise as RMS. By default, the data system uses peak-to-peak for the noise calculation. RMS is automatically selected if you determine the noise region manually. |
| Min Peak Width | Type the minimum number of scans required in a peak. The valid range is 0 to 100 scans. The default value is 3 scans. This value is used by the ICIS peak detection algorithm. |
| Multiplet Resolution | Type the minimum separation in scans between the apexes of two potential peaks. This is a criteria to determine if two peaks are resolved. The valid range is 1 to 500 scans. The default value is 10 scans. This value is used by the ICIS peak detection algorithm. |

Table 43. ICIS Advanced page parameters (Sheet 2 of 2)

| Parameter | Description |
|---------------------|---|
| Area Tail Extension | Type the number of scans past the peak endpoint to use in averaging the intensity. The valid range is 0 to 100 scans. The default value is 5 scans. This value is used by the ICIS peak detection algorithm. |
| Area Scan Window | Type the number of allowable scans on each side of the peak apex. The valid range is 0 to 100 scans. The default value of 0 scans specifies that all scans from peak start to peak end are to be included in the area integration. This value is used by the ICIS peak detection algorithm. |

Flags Page – User Identification Settings Dialog Box

The Flags page of the User Identification Settings dialog box shows the current detection flagging thresholds in use for the selected compound. These values are used in determining if the peak detection is within user-specified limits. They do not alter the way calculations are made. An entered value of 0.0 forces the flag to be false.

This table describes the parameters on the Flags page of the User Identification Settings dialog box.

Table 44. Flags page parameters – User Identification Settings dialog box

| Parameter | Description |
|------------------|---|
| Area Threshold | View or set a value for the current Area Threshold (AT) flag. The data system sets the AT flag in the result file if the quantified peak has an area that is lower than the entered value. |
| Height Threshold | View or set a value for the current Height Threshold (HT) flag. The data system sets the HT flag in the result file if the quantified peak has a height that is lower than the entered value. |

View Sample Types Dialog Box

When opening Quan Browser, the Xcalibur data system first checks to confirm that the file you select is valid. After verifying that all the files exist and can be opened, but before displaying any data, the View Sample Types dialog box opens to prompt you to display only “Standards and QC samples” or “All samples.”

The Don't Ask Again check box is provided so that you do not have to see the dialog box again after making your initial selection.

- If you select the Don't Ask Again check box, choose **Options > Enable Warnings** from the Quan Browser menu bar to reset the display of this dialog box and all other message type dialog boxes.
- If you do not select the Don't Ask Again check box, the Enable Warnings menu command is not available.

This table describes the parameters in the View Sample Types dialog box.

Table 45. View Sample Types dialog box parameters

| Parameter | Description |
|------------------------|---|
| Viewing Options | |
| Show Standards and QCs | Display only Standards and QCs in the Quan Browser results grid view. The data system does not display blanks and Unknowns. Select one of these tabs: Standards or QCs. |
| Show All Sample Types | Display Standards, QCs, Blanks, and Unknowns in the Quan Browser grid view. You can select from the following tabs: All, Standards, QCs, Blanks, or Unknowns. |
| Don't Ask Again | Decide whether you want to see the current message box or dialog box in the future. For example, if you always select the Show All Sample Types option and never select the Show Standards or QCs option, you might want to turn off the View Sample Types dialog box. When you select the Don't Ask Again check box, the Quan Browser application does not display the View Samples dialog box each time you open a new sequence. However, after you close the Quan Browser application, the data system automatically clears the Don't Ask Again check box, and the next time you open the Quan Browser application, the View Sample Types box opens. To turn on the display of all message boxes and dialog boxes, choose Options > Enable Warnings . |

Index

A

accessing the manuals as PDF files [viii](#)
advanced integration parameters [43](#), [129](#)
Avalon Integration page [142](#)

B

blue baselines [46](#)
bracket types
 Non-overlapping [14](#)
 Open [14](#)
 Overlapping [14](#)
 Quan Browser overview [12](#)
 unbracketed [12](#)
Brackets/Groups In Use box [18](#)

C

calibration curve
 modifying parameters [54](#)
 replicates [11](#)
 using external standard (figure) [7](#)
calibration file, using [12](#)
Calibration Settings dialog box [98](#)
chromatogram plot view
 description [19](#)
 reviewing [33](#)
 working in [33](#)
chromatograms
 integrating peaks automatically [41](#)
 integrating peaks manually [46](#)
Columns command [31](#)
component list, setting [17](#)
contacting us [ix](#)
continuing calibration method [12](#)

D

data points, excluding from the calibration curve [59](#)
detection limit [5](#)

Detection page, User Identification Settings dialog box [40](#)
documentation, Xcalibur data system [vii](#)

E

excluding data points [59](#)
external calibration file [12](#)
external standards, considering variables for [6–7](#)

F

File menu [76](#)
file types supported [15](#), [50](#), [124–125](#)

G

Genesis Advanced page [43](#), [145](#)
Genesis Integration page [139](#)
Go To menu [80](#)

H

Help menu [80](#)

I

ICIS Advanced page [147](#)
ICIS Integration page [141](#)
Identification page, User Identification Settings dialog box [38](#)
integration
 manual [46](#)
 parameters advanced [67](#)
integration algorithms
 Avalon [142](#)
 Genesis [139](#), [145](#)
 ICIS [141](#), [147](#)
internal standards (ISTDs)
 choosing [9](#)
 considering variables for [8](#)
 definition [8](#)
 using, for quantitation [8](#)

K

Keys box, comments in 40

L

Levels page, Calibration Settings dialog box 57

limit

of detection 5

of quantitation 5

lower quantitation limit 5

M

manual peak integration 46

N

non-overlapped, bracket type 14

O

open, bracket type 14

Options menu 79

overlapped, bracket type 14

P

Peak Detection Settings command, User Identification Settings dialog box 37

Peak Information dialog box 35

peaks, chromatographic

integrating automatically 46

integrating manually in chromatogram 46

processing methods, saving 23

Q

Qual Browser window, sending result file to 48

Quan Browser dialog boxes

Calibration Settings dialog box

Curve page 98

Flags page 100

Isotope% page 101

Levels page 104

list of pages 98

Type page 105

Peak Information dialog box

Chro page 115

description 108

Flags page 111

Info page 110

Info/More Info page 116

list of pages 108

More Flags page 114

More Info page 117

No Peak page 109

opening 35

Spectrum page 118

Suitability page 118

Quantitation Results Sorting Order dialog box 119

Reports dialog box 68, 121

Result List Column Hiding dialog box 31, 126

Select Level dialog box 127

Select Report Samples dialog box 128

User Identification Settings dialog box 37, 129

View Sample Types dialog box 15

Quan Browser views

calibration curve view 92

chromatogram plot view 88

components list 88

results grid 84

spectrum plot view 90

Quan Browser window

chromatogram view 19

components list 17

elements 74

features 17

menu bar 76

opening 15

opening files 15

Quantitation Results Sorting Order dialog box 32

title bar 75

quanbrowser files 21

quantitation limits 5

quantitation range 5

Quantitation Results Sorting Order dialog box 32, 119

quantitative analysis

definition 2

sources of error 8

using internal standards for 8

R

Replace Calibration command 18

replicates 11

reports

file types supported 124–125

generating in Quan Browser 68

Reports dialog box

opening 68

parameter descriptions 121

Reset Scaling command 49

result files, saving 21

Result List Column Hiding dialog box 31, 126

results grid
 changing the sort order 32
 column descriptions 85
 column headings 31
 displaying columns 31
 editing a sequence 25
 hiding columns 31
 shortcut menu 87
 working in 18

S

saving
 changes in Quan Browser 26
 QuanBrowser files 21
 user identification settings 38
 scaling, resetting 49
 Select Level dialog box 127
 Select Report Samples dialog box 128
 sending a result file to Qual Browser 48
 Set Sorting Order command 32
 Show Calibration Curve command 49
 Show Peak Info command 35
 Show Standards and QC commands 15
 sort order 32
 Spectrum at Peak Apex command 49
 Spectrum at Peak Left Edge command 49
 Spectrum at Peak Right Edge command 49
 Spectrum Plot view 48
 spreadsheets, exporting the results grid to 23
 standard
 clear 12
 update 12
 system suitability results 36

T

target component 5
 toolbar, Quan Browser 81

U

unbracketed sequence 12
 upper quantitation limit 5
 user guides vii

User Identification Settings dialog box
 description 129
 Detection page 40
 ICIS Advanced page 43
 ICIS Integration page 41
 Identification page 38, 130
 User Peak Detection Settings command 37

V

variables, discussion of
 quantitation with external standards 6
 quantitation with internal standards 8
 View menu 78
 View Sample Types dialog box 15
 View Spectrum Plot command 48

W

Warning dialog box 15
 working in the results grid 18

X

Xcal files 12
 Xcalibur data system
 acquiring and processing data with, overview 2
 QuanBrowser file, saving 21
 quantitative analysis procedure 2

Z

Zoom menu, Quan Browser 79

