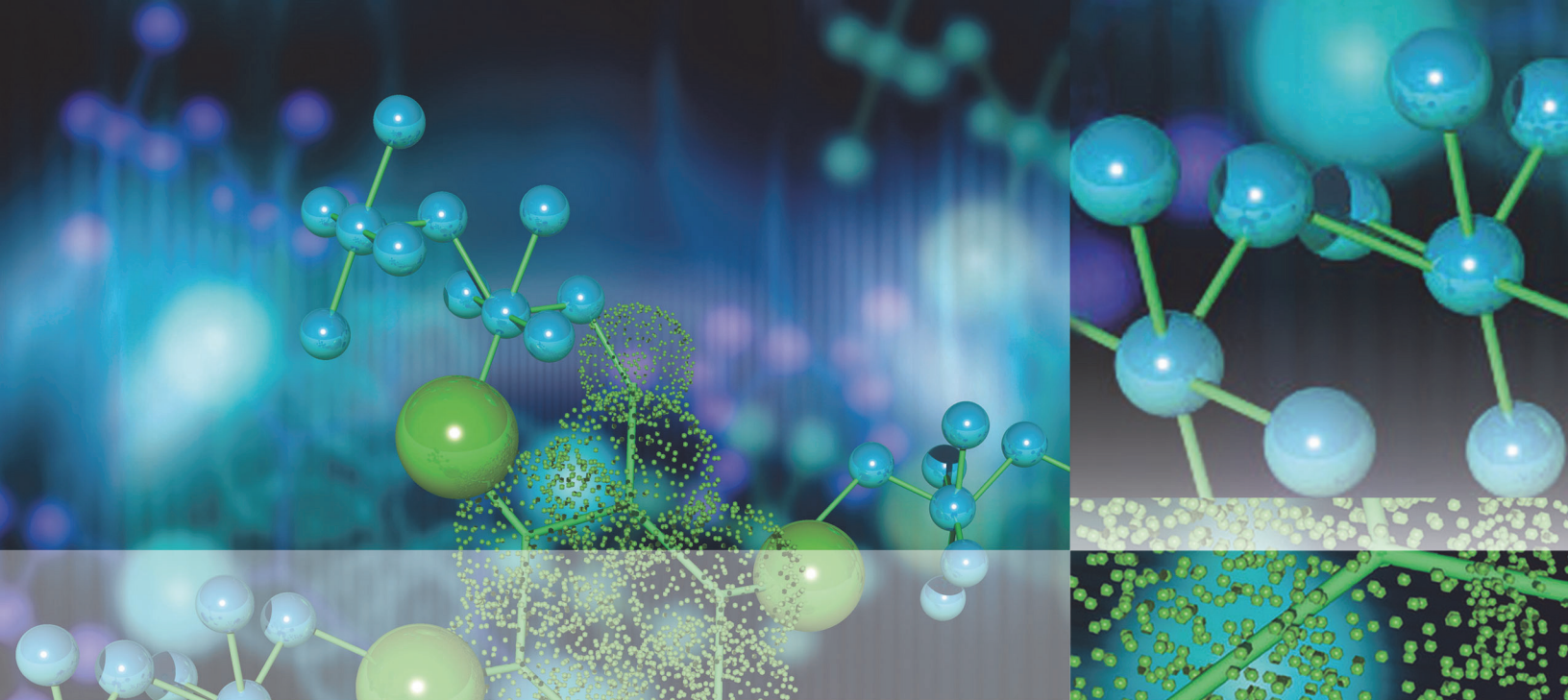




Xcalibur

LCquan

Quantitative Analysis of a Three-Drugs Data Set

Tutorial

Software Version 3.0

XCALI-97816 Revision A April 2016

Thermo
SCIENTIFIC

© 2016 Thermo Fisher Scientific Inc. All rights reserved.

DCMS^{Link}, Foundation, and LCquan are trademarks, and Exactive, Thermo Scientific, TSQ Endura, TSQ Quantiva, and Xcalibur are registered trademarks of Thermo Fisher Scientific Inc. in the United States.

Excel, Microsoft, and Windows are registered trademarks of Microsoft Corporation 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, April 2016

Software version: (Thermo) Foundation 3.1.100, Xcalibur 4.0, LC Devices 3.0 QF1 or later, DCMS^{Link} 2.14, (Microsoft) Windows 7 Professional SP1 (32-bit or 64-bit) and Office 2010, one of the following mass spectrometer (MS) drivers: TSQ Endura[™] 2.0 and 2.1, TSQ Quantiva[™] 2.0 and 2.1, Exactive Series 2.7

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

Contents

	Preface	v
	Accessing Documentation	v
	Special Notices	vi
	Contacting Us	vi
Chapter 1	Tutorial Exercises	1
	Introduction	1
	Exercise 1: Creating a New Study and Workbook	2
	Exercise 2: Creating an Acquisition Sequence	7
	Using the New Acquisition Sequence Wizard	8
	Editing the Acquisition Sequence	14
	Naming Components and Assigning Calibration Levels	17
	Exercise 3: Exploring the Data	19
	Opening a Raw Data File and Displaying the Chromatogram	20
	Exploring Peak Detection and Integration Parameters	21
	Reviewing the Explore Results and Determining Major Product Ions	23
	Exercise 4: Creating a Processing Method	27
	Opening the Create Method Page and Using the New Method Wizard	27
	Identifying the Internal Standard	29
	Specifying Calibration Curve Parameters for the Target Compounds	30
	Entering Component Identification Parameters and Integrating the Peaks	32
	Saving the Processing Method	42
	Saving the Workbook and Exiting the LCquan Application	42
	Exercise 5: Creating a Processing Sequence	42
	Opening the Create Sequence Page and Using the New Processing Sequence Wizard	42
	Associating Raw Data Files with Sample Types	43
	Reviewing the Processing Sequence	48
	Exercise 6: Processing the Raw Data Files and Reviewing the Analytical Results	49
	Processing the Raw Data Files	50
	Reviewing the Calibration Standards	51
	Reviewing the Unknowns	54
	Exercise 7: Generating a Custom Report	57
	Saving the Workbook and Exiting the LCquan Application	59
	Index	61

Preface

The LCquan™ application is part of the Thermo Xcalibur™ mass spectrometry data system. This tutorial provides step-by-step procedures to perform quantitative analysis with sample data.

Contents

- [Accessing Documentation](#)
- [Special Notices](#)
- [Contacting Us](#)

❖ To suggest changes to the documentation or to the Help

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



Accessing Documentation

The LCquan application includes complete documentation. In addition to this guide, you can also access the following documents as PDF files from the data system computer:

- *Foundation Administrator Guide*
- *LCquan User Guide*

❖ To view the product manuals

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

❖ **To view user documentation from the Thermo Fisher Scientific website**

1. Go to thermofisher.com.
2. Click the **Services & Support** tab.
3. On the right, click **Manuals & Protocols**.
4. In the Refine Your Search box, search by the product name.
5. From the results list, click the title to open the document in your web browser, save it, or print it.

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

Special Notices

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

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

Note Highlights information of general interest.

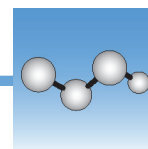
Tip Highlights helpful information that can make a task easier.

Contacting Us

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

Contact us	Customer Service and Sales	Technical Support
	(U.S.) 1 (800) 532-4752	(U.S.) 1 (800) 532-4752
	(U.S.) 1 (561) 688-8731	(U.S.) 1 (561) 688-8736

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



Tutorial Exercises

This tutorial contains a step-by-step procedure for performing quantitative analysis on the example data set provided in the Xcalibur\QuanRoot\Example Study folder of your Xcalibur data system.

Contents

- [Introduction](#)
- [Exercise 1: Creating a New Study and Workbook](#)
- [Exercise 2: Creating an Acquisition Sequence](#)
- [Exercise 3: Exploring the Data](#)
- [Exercise 4: Creating a Processing Method](#)
- [Exercise 5: Creating a Processing Sequence](#)
- [Exercise 6: Processing the Raw Data Files and Reviewing the Analytical Results](#)
- [Exercise 7: Generating a Custom Report](#)
- [Saving the Workbook and Exiting the LCquan Application](#)

Introduction

The data used in this tutorial comes from the analysis performed on a mixture of drugs in the applications laboratories at Thermo Fisher Scientific by using LC/MS/MS techniques in the electrospray mode.

The three-drugs mixture contains paroxetine, nefazodone, and alprazolam as the target compounds. In the example, deuterated alprazolam (alprazolam-d5) functions as an internal standard (ISTD). The internal standard in the calibration standards has a concentration of 1.0 ng/ml. Calibration standards have levels with amounts of 25, 50, 75, 100, 250, 500, 750, and 1000 pg/ml.

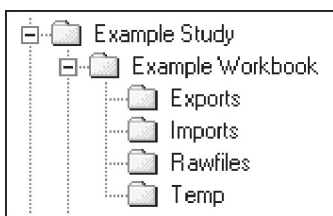
Note

- To open a completed three-drugs study and workbook, choose **Open an Existing Workbook** on the LCQuan startup page, browse to ...\\Xcalibur\\QuanRoot\\ Example Study\\3-drugs Workbook, and click **OK**.
- For a more complete description of LCQuan features, refer to the *LCQuan User Guide*.

Exercise 1: Creating a New Study and Workbook

When you start a new quantitative analysis project, the LCQuan application creates a hierarchical folder structure for the project on the hard drive, as shown in [Figure 1](#).

Figure 1. Folder structure for the LCQuan application



Use the Study folder, directly beneath the quantitation root in the file structure, to organize one or more workbook folders. Each workbook folder corresponds to an individual quantitative analysis project. The workbook folder contains the LCQuan file (.lqn), the Instrument Method files (.meth), and the audit log files (.mdb). In addition, the workbook folder contains the following folders: Exports, Imports, Rawfiles, and Temp.

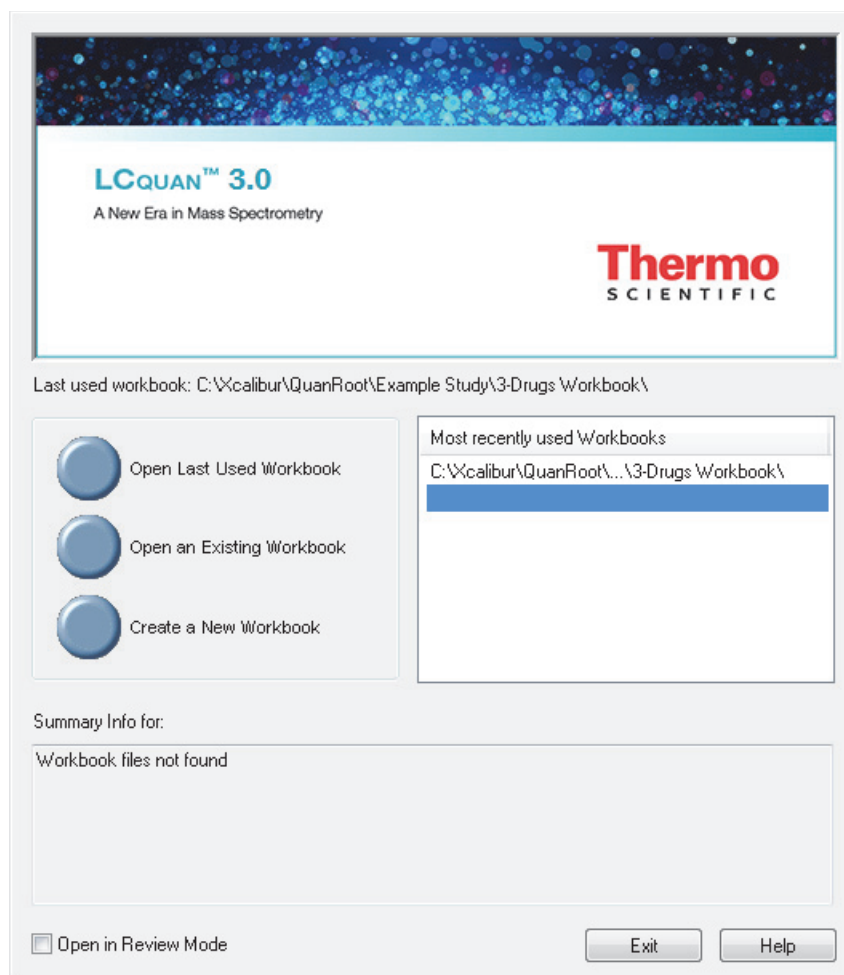
To start the quantitative analysis of the three-drugs data, create a new study and workbook before importing raw data files.

❖ To create the new study and workbook

1. Choose **Start > All Programs > Thermo Xcalibur > LCQuan** to open the LCQuan application (see [Figure 2](#)).

From the LCQuan startup dialog box, you can open the last used workbook, open an existing workbook, or create a new workbook.

Figure 2. LCquan open workbook window



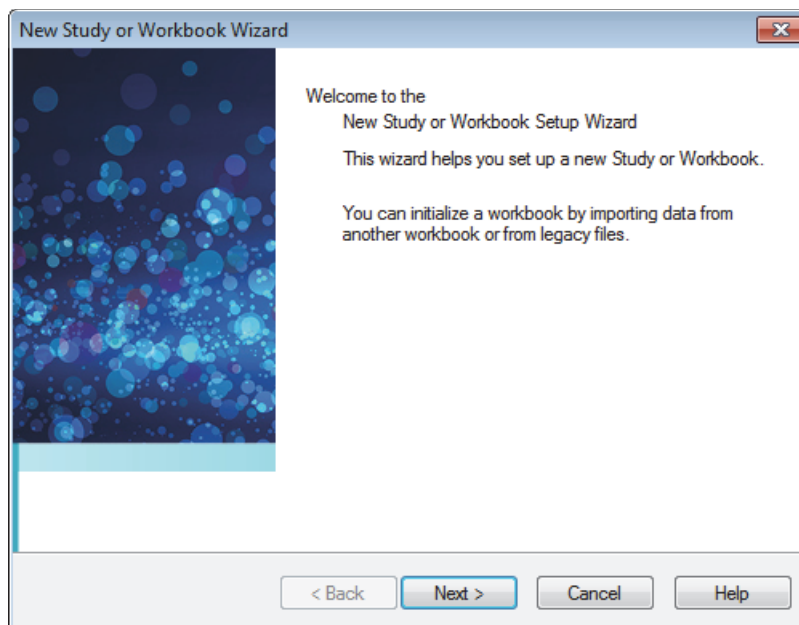
2. Ensure that the **Open in Review Mode** check box remains cleared (default).
3. Click **Create a New Workbook**.

The welcome page of the New Study or Workbook Wizard opens (see [Figure 3](#)).

1 Tutorial Exercises

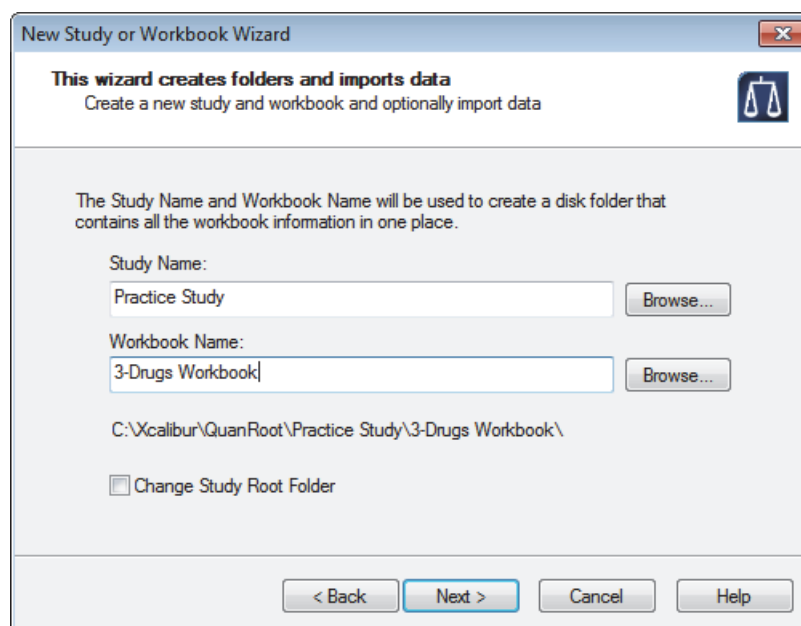
Exercise 1: Creating a New Study and Workbook

Figure 3. Welcome page of the New Study or Workbook Wizard



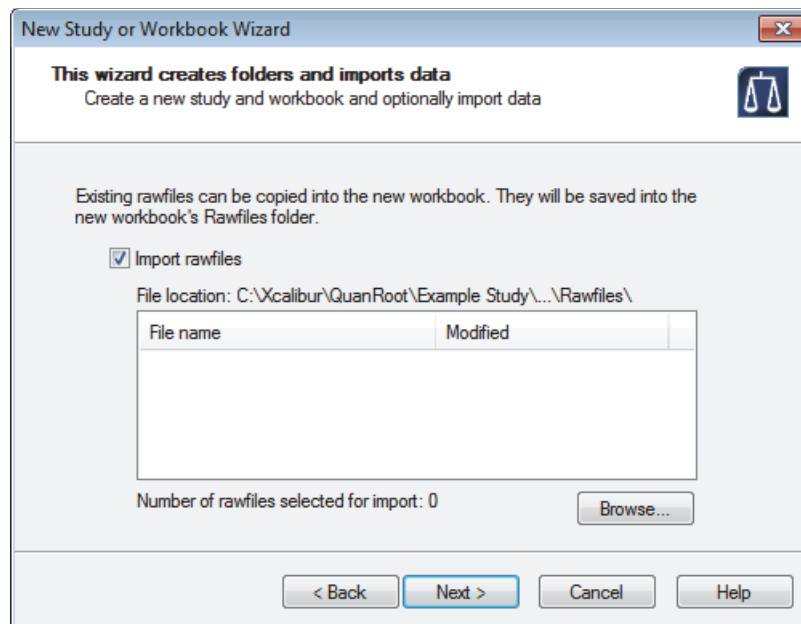
4. Click **Next**, and then name the study and workbook folders, as follows (see [Figure 4](#)):
 - a. In the Study Name box, type **Practice Study**.
 - b. In the Workbook Name box, type **3-Drugs Workbook**.

Figure 4. Create new study and workbook page



5. Click **Next**, and then select the raw data files to import into the new workbook, as follows (see [Figure 5](#)):
 - a. Select the **Import Rawfiles** check box.
 - b. Click **Browse**.

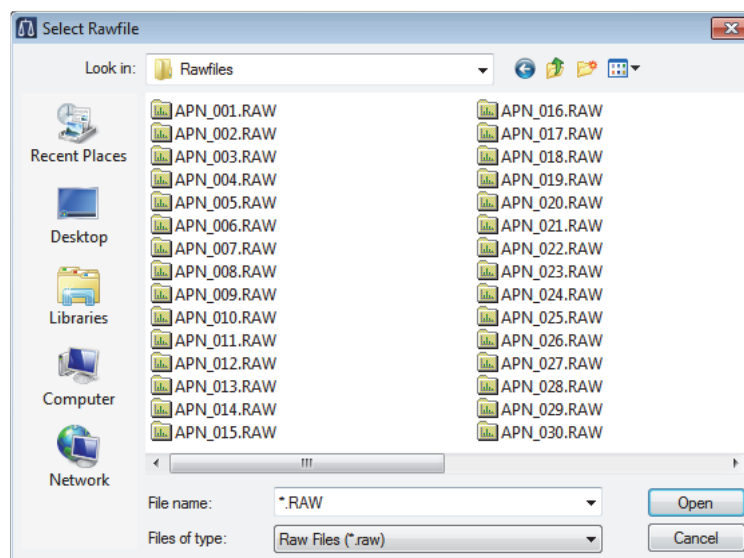
Figure 5. Import raw data files page



- c. In the Select Rawfile dialog box, browse to the Rawfiles folder:
... \Xcalibur\QuanRoot\Example Study\3-Drugs Workbook\Rawfiles

[Figure 6](#) shows the content of the Rawfiles folder.

Figure 6. Select Rawfile dialog box



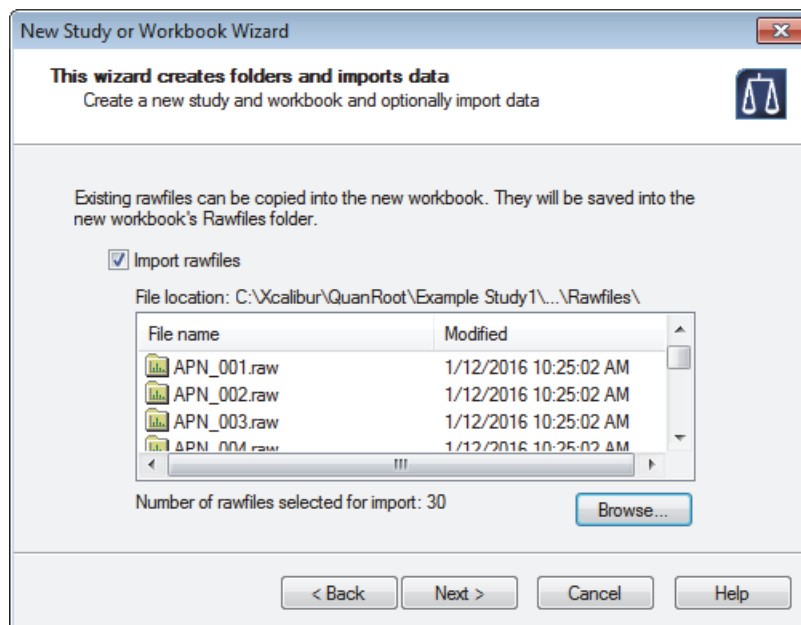
1 Tutorial Exercises

Exercise 1: Creating a New Study and Workbook

- d. To select all 59 raw data files, select **APN_001**, press SHIFT, and select **APN_59**.
- e. Click **Open** to accept your selection.

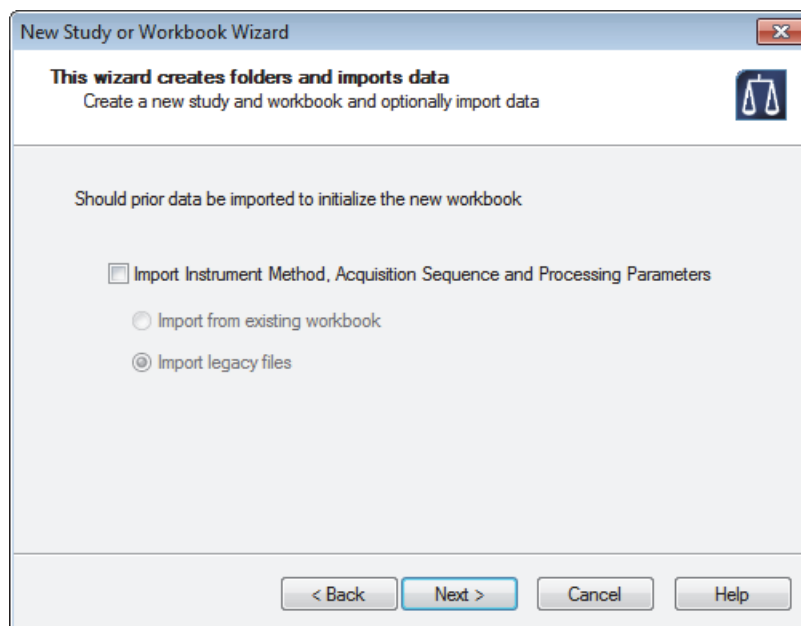
The wizard displays the selected raw data files, as shown in [Figure 7](#).

Figure 7. Import rawfiles page



6. Click **Next**.
7. Clear the **Import Instrument Method, Acquisition Sequence and Processing Parameters** check box (see [Figure 8](#)).
8. Select either the **Import from Existing Workbook** or the **Import Legacy Files** option.

Figure 8. Import data page



9. Click **Next**.
10. On the final New Study or Workbook Wizard page, click **Finish**.

The LCquan application creates the new study and workbook folders and imports the raw data files into the Rawfiles folder.

After you create a new workbook, the LCquan application opens the Instrument Setup view. Because the data already exists, go to [“Exercise 2: Creating an Acquisition Sequence.”](#)

Saving the Workbook and Exiting the LCquan Application

If you want to take a break, you can save your workbook and open it again when you are ready to continue. For instructions about saving a workbook, see [“Saving the Workbook and Exiting the LCquan Application”](#) on page 59.

Exercise 2: Creating an Acquisition Sequence

The acquisition sequence contains the list of samples and defined settings that the LCquan application uses to acquire data. The acquisition sequence can contain unknown samples, calibration standard samples, quality control samples, or blank samples. Use the Acquisition view to define an acquisition sequence, run individual samples and sequences, and monitor real-time data acquisition.

This exercise uses a wizard to define an acquisition sequence for the quantitative analysis of the three-drugs data.

Using the New Acquisition Sequence Wizard

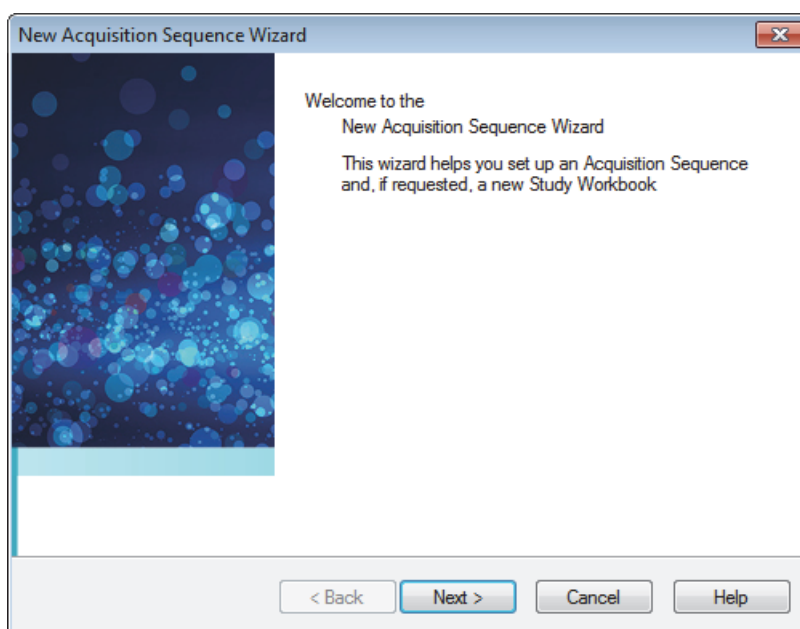
❖ To define an acquisition sequence for the three-drugs example

1. In the navigation pane, click .

Because no acquisition sequence is defined in the workbook, the New Acquisition Sequence Wizard opens (see [Figure 9](#)).

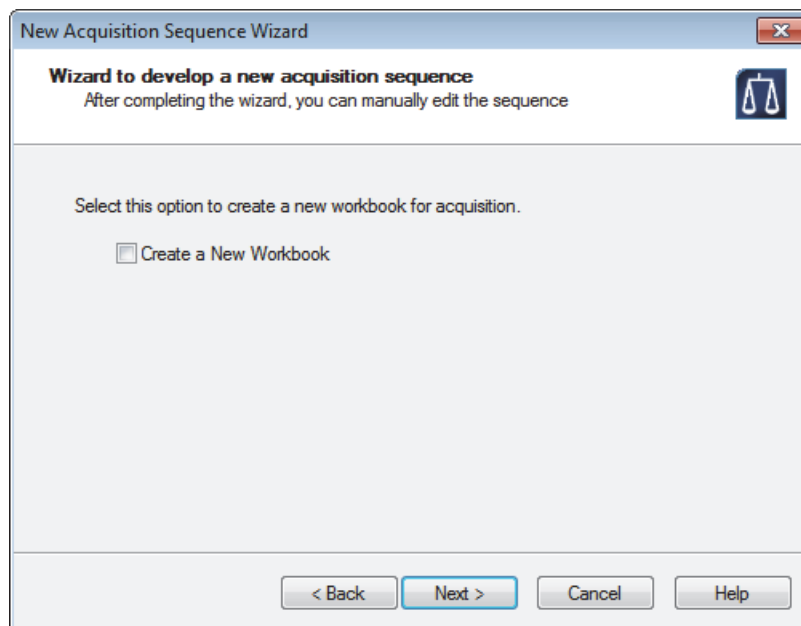
Tip If the New Acquisition Sequence Wizard does not appear, right-click the sequence grid and choose **New Acquisition Sequence Wizard**.

Figure 9. Welcome page of the New Acquisition Sequence Wizard



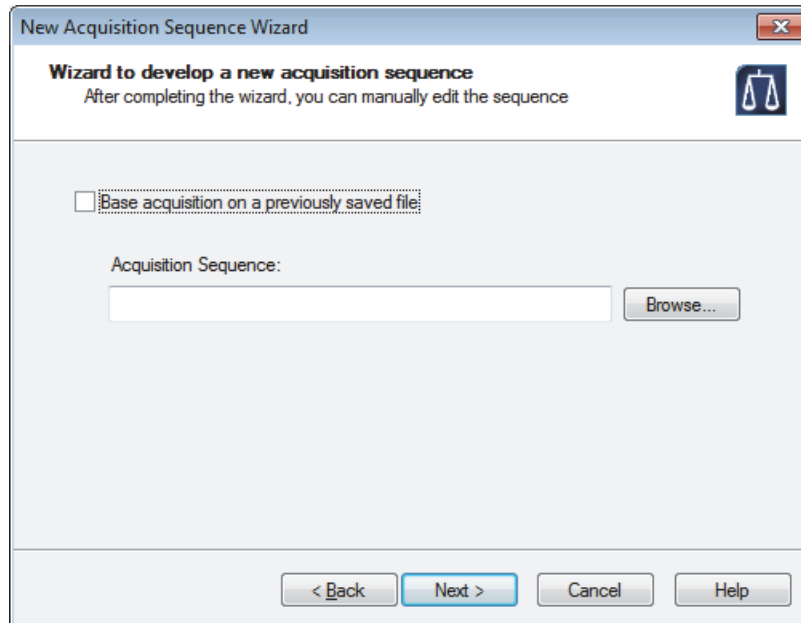
2. Click **Next**, and on the next page ensure that the **Create a New Workbook** check box is cleared (see [Figure 10](#)).

Figure 10. Create a new workbook page



3. Click **Next**, and on the next page ensure that the **Base Acquisition on a Previously Saved File** check box is cleared (see [Figure 11](#)).

Figure 11. Acquisition sequence page



1 Tutorial Exercises

Exercise 2: Creating an Acquisition Sequence

4. Click **Next**, and then specify the raw data file names (see [Figure 12](#)):

- Base file name: **APN_**
- Starting number: **1**

Figure 12. Sample sequence name page

New Acquisition Sequence Wizard

Wizard to develop a new acquisition sequence
After completing the wizard, you can manually edit the sequence

Names are generated for the samples by using a base name and appending a numerical value that is incremented for each sample. The base name and starting numerical value are entered here.

An example might be: Base file name = "TestSample", Starting number 100. The first file name would be TestSample100, the next file name would be TestSample101 ...

Base file name:

Starting number:

< Back Next > Cancel Help

5. Click **Next**.

The LCQuan application appends a three-digit number (001-999) to the base file name so that each sample in a sequence has a unique identification.

6. Specify the unknown sample information exactly as indicated in [Figure 13](#).

Figure 13. Sample information page

The screenshot shows the 'New Acquisition Sequence Wizard' dialog box. The title bar reads 'New Acquisition Sequence Wizard'. Below the title bar, the text says 'Wizard to develop a new acquisition sequence' and 'After completing the wizard, you can manually edit the sequence'. The main area contains the instruction 'Specify how many samples to create and what they will be named.' followed by three input fields: 'Number of Samples:' with the value '4', 'Injections per Sample:' with the value '1', and 'Base Sample ID:' with the value 'APN_'. At the bottom, there are four buttons: '< Back', 'Next >', 'Cancel', and 'Help'.

7. Click **Next** (see [Figure 14](#)).

The LCQuan application generates sample IDs that uniquely identify the unknown samples in the sequence by appending numerical values to a base sample ID.

Figure 14. Tray and vial information page

The screenshot shows the 'New Acquisition Sequence Wizard' dialog box, specifically the 'Tray and vial information' page. The title bar reads 'New Acquisition Sequence Wizard'. Below the title bar, the text says 'Wizard to develop a new acquisition sequence' and 'After completing the wizard, you can manually edit the sequence'. The main area contains the instruction 'Specify the vial positions for each of the samples based on the type of autosampler tray in use.' followed by four input fields: 'Tray Type:' with a dropdown menu, 'Initial Vial Position:' with the value '1', 'Re-Use Vial Positions' with a checked checkbox, and 'Injection Volume obtained from Autosampler Method' with a checked checkbox. Below these, there is an 'Injection Volume' field with the value '10.00'. At the bottom, there are four buttons: '< Back', 'Next >', 'Cancel', and 'Help'.

8. Leave the autosampler settings as they are.

1 Tutorial Exercises

Exercise 2: Creating an Acquisition Sequence

9. Click **Next**, and then specify the calibration standards exactly as indicated in [Figure 15](#).

Figure 15. Add standards page

The screenshot shows the 'New Acquisition Sequence Wizard' dialog box. The title bar reads 'New Acquisition Sequence Wizard'. Below the title bar, the text says 'Wizard to develop a new acquisition sequence' and 'After completing the wizard, you can manually edit the sequence'. There is a balance scale icon in the top right corner. The main area contains the following options:

- ☒ Add Standards
 - ☐ Based on existing cal levels from method
 - ☒ Based on automatically generated cal levels
 - Number of cal levels: 8
 - Cal level base name: Cal
- Number of Injections per Level: 10
- ☒ Add Blanks
- ☒ Fill in Sample ID for Standards

At the bottom, there are four buttons: '< Back', 'Next >', 'Cancel', and 'Help'.

10. Click **Next**, and then specify the quality control options exactly as indicated in [Figure 16](#).

The LCQuan application generates names for the levels by appending numerical values to the base name.

Figure 16. Add QCs page

The screenshot shows the 'New Acquisition Sequence Wizard' dialog box. The title bar reads 'New Acquisition Sequence Wizard'. Below the title bar, the text says 'Wizard to develop a new acquisition sequence' and 'After completing the wizard, you can manually edit the sequence'. There is a balance scale icon in the top right corner. The main area contains the following options:

- ☒ Add QCs
 - ☐ Based on existing QC levels from method
 - ☒ Based on automatically generated QC levels
 - Number of QC levels: 3
 - QC level base name: QC_
- ☒ Add Blanks
- ☒ Fill in Sample ID for QCs

At the bottom, there are four buttons: '< Back', 'Next >', 'Cancel', and 'Help'.

11. Click **Next**, and then specify temporary component names for the target drugs exactly as shown in [Figure 17](#).

Until you name your target components, the system temporarily names the components *Component1-4*.

Figure 17. Component names page

The screenshot shows the 'New Acquisition Sequence Wizard' dialog box. The title bar reads 'New Acquisition Sequence Wizard'. The main heading is 'Wizard to develop a new acquisition sequence' with a subtitle 'After completing the wizard, you can manually edit the sequence'. There is a checkbox labeled 'Use processing method component names' which is currently unchecked. Below this, there is a text input field for 'Number of components' containing the value '4', and another text input field for 'Component base name' containing the text 'Component'. At the bottom of the dialog are four buttons: '< Back', 'Next >', 'Cancel', and 'Help'.

12. Click **Next**, and then accept the default user labels and values (see [Figure 18](#)).

These labels and values are used in the header of the raw data file.

Figure 18. User labels and values page

The screenshot shows the 'New Acquisition Sequence Wizard' dialog box, specifically the 'User labels and values' page. The title bar and main heading are the same as in Figure 17. Below the heading, there is a paragraph: 'When creating a new sequence, user configurable labels can be applied to each sample. The following will be used when creating new samples in the sequence.' Below this paragraph are five rows of input fields. Each row has a 'User Label' text input and a 'User Value' dropdown menu. The labels are: 'User Label 1: Study', 'User Label 2: Client', 'User Label 3: Laboratory', 'User Label 4: Company', and 'User Label 5: Phone'. The corresponding values in the dropdowns are: 'User Value 1: \$Study', and the others are empty. At the bottom are the same four buttons: '< Back', 'Next >', 'Cancel', and 'Help'.

13. Click **Next** and then click **Finish** to close the wizard.

The Acquisition – Setup Sequence page opens (see [Figure 19](#)).

1 Tutorial Exercises

Exercise 2: Creating an Acquisition Sequence

Figure 19. Acquisition – Setup Sequence page

	Sample Type	Filename	Sample ID	Level	ISTD Coef	Dil Factor	Vial Pos	Inj Vol	Sample	S
1	Blank	APN_001		NA	0.000	1.000	A:1	From AS	0.000	0.1
2	Standard	APN_002	APN_002	Cal_01	0.000	1.000	1	From AS	0.000	0.1
3	Standard	APN_003	APN_003	Cal_01	0.000	1.000	1	From AS	0.000	0.1
4	Standard	APN_004	APN_004	Cal_01	0.000	1.000	1	From AS	0.000	0.1
5	Standard	APN_005	APN_005	Cal_01	0.000	1.000	1	From AS	0.000	0.1
6	Standard	APN_006	APN_006	Cal_01	0.000	1.000	1	From AS	0.000	0.1
7	Standard	APN_007	APN_007	Cal_02	0.000	1.000	2	From AS	0.000	0.1
8	Standard	APN_008	APN_008	Cal_02	0.000	1.000	2	From AS	0.000	0.1
9	Standard	APN_009	APN_009	Cal_02	0.000	1.000	2	From AS	0.000	0.1
10	Standard	APN_010	APN_010	Cal_02	0.000	1.000	2	From AS	0.000	0.1
11	Standard	APN_011	APN_011	Cal_02	0.000	1.000	2	From AS	0.000	0.1
12	Standard	APN_012	APN_012	Cal_03	0.000	1.000	3	From AS	0.000	0.1
13	Standard	APN_013	APN_013	Cal_03	0.000	1.000	3	From AS	0.000	0.1
14	Standard	APN_014	APN_014	Cal_03	0.000	1.000	3	From AS	0.000	0.1
15	Standard	APN_015	APN_015	Cal_03	0.000	1.000	3	From AS	0.000	0.1
16	Standard	APN_016	APN_016	Cal_03	0.000	1.000	3	From AS	0.000	0.1
17	Standard	APN_017	APN_017	Cal_04	0.000	1.000	4	From AS	0.000	0.1
18	Standard	APN_018	APN_018	Cal_04	0.000	1.000	4	From AS	0.000	0.1
19	Standard	APN_019	APN_019	Cal_04	0.000	1.000	4	From AS	0.000	0.1
20	Standard	APN_020	APN_020	Cal_04	0.000	1.000	4	From AS	0.000	0.1
21	Standard	APN_021	APN_021	Cal_04	0.000	1.000	4	From AS	0.000	0.1
22	Standard	APN_022	APN_022	Cal_05	0.000	1.000	5	From AS	0.000	0.1

Editing the Acquisition Sequence

Although you do not acquire data in this example, you can edit the acquisition sequence as if you were. This acquisition sequence produces the raw data files in the data set. You can use this acquisition sequence as the processing sequence later on.

❖ To edit the acquisition sequence

1. Delete rows 51 through 92, as follows:
 - a. Select the **51** in the number column at the left of the row and drag straight down to row **92**.
 - b. Right-click the highlighted rows and choose **Delete Rows**.

The selected rows are deleted.

2. Insert additional empty rows before each of the rows with the Blank sample type 1, 42, and 46, as follows:

- a. Click the first blank row (it turns turquoise), right-click, and choose **Insert Rows**.

An empty row appears above the selected row (see Figure 20).

Repeat this until you have three new empty rows above the original Blank sample type row.

Figure 20. New empty rows

	Sample Type	Filename	Sample ID
1			
2			
3			
4	Blank	APN_001	

- b. To copy information into these empty rows, click the number of the original Blank row to select the row, right-click, and choose **Duplicate Selected Samples**.

1			
2			
3			
4	Blank	APN_001	

Selected row

Note You can select a row only by clicking its number in the far left column.

The duplicated row is copied to the end of the sequence, and the sequence display scrolls down to this added row.

- c. To move the information in the added row (at the end of the sequence) into one of your empty rows (at the beginning of the sequence), drag the number of the added row to the top of the sequence grid and drop the row.

The row you moved overwrites the row **below** the red bar (see Figure 21).

Figure 21. Placement indicator of moved row

	Sample Type	Filename	Sample ID
1			
2			
3			
4	Blank	APN_001	APN001

Repeat steps b and c until you have moved the information into the three new empty rows (see Figure 22). Do not worry about the duplicate file names. You will fix those later.

Figure 22. Copied blank rows

	Sample Type	Filename	Sample ID
1	Blank	APN_001	
2	Blank	APN_001	
3	Blank	APN_001	
4	Blank	APN_001	

1 Tutorial Exercises

Exercise 2: Creating an Acquisition Sequence

- d. Repeat steps **a** through **c** for each of the original rows: 1, 42, and 46.

Note Now that you have added rows above the original rows, the numbers have changed. Make sure you copy the Blank rows that were originally numbers 42 and 46.

3. For the newly added rows, change the Sample Type to QC, as follows:

- a. In each of the new rows, click the Sample Type column.

A down arrow appears.

- b. From the list, select **QC**.

The remaining quality control information automatically populates the row.

Figure 23. Quality control sample type

	Sample Type	Filename	Sample ID	Level
1	QC	APN_001		QC_01
2	Standard	APN_001		QC_01
3	QC	APN_001		QC_01
4	Blank	APN_001		NA
5	Unknown Standard	APN_002	APN_002	Cal01

- c. Repeat steps **a** and **b** for the remaining new rows.

Remember that you created a total of nine new rows.

4. Change the Levels to QC_*, as follows:

- a. Change the first three new QC row Levels to **QC_01**.
- b. Change the next three new QC row Levels to **QC_02**.
- c. Change the last three new QC row Levels to **QC_03**.

5. Update the file names, as follows:
 - a. Right-click the FileName column and choose **Fill Down** (see Figure 24).

Figure 24. Fill Down dialog box

Select Columns		
☐ Sample Type	☐ Sample Name	☐ Aria LC Loop Fill
☒ FileName	☐ Study	☐ Aria LC Flow Rate
☐ Sample ID	☐ Client	☐ Aria LX/TX
☐ Level	☐ Laboratory	
☐ ISTD Corr Amt	☐ Company Name	
☐ Dil Factor	☐ Phone	
☐ Vial Pos	☐ Comment	
☐ Inj Vol	☐ Barcode	
☐ Sample Vol	☐ Aria MCM1	
☐ Sample Vt	☐ Aria MCM2	

Fill From Row To Row Using Row

☐ Create samples in selected empty rows during fill down

OK Select All Clear All Cancel Help

- b. In the Fill Down dialog box and with the Filename column selected, specify the rows you want to fill:

Fill From Row **1** To Row **59**

- c. Specify the number of the row you want to duplicate:
 - Using Row **1**
 - d. Clear the **Create Samples in Selected Empty Rows During Fill Down** check box.
 - e. Click **OK**.

The system rennumbers the filenames in consecutive order.

6. Choose **File > Save** to save the acquisition sequence and workbook.

Naming Components and Assigning Calibration Levels

Use the Acquisition Levels dialog box to change the component names and define or modify the calibration levels and QC levels used by the acquisition sequence.

❖ To name components and assign calibration levels

1. Right-click the sequence grid on the Setup Sequence page and choose **Standard and QC Levels**.

The Acquisition Levels dialog box opens (see [Figure 25](#)).

2. Change the component names by selecting the default name in the Components table and entering the new name:
 - a. Change Component1 to **Paroxetine**.
 - b. Change Component2 to **Alprazolam**.
 - c. Change Component3 to **Alprazolam-d5**.
 - d. Change Component4 to **Nefazodone**.
 - e. To save the new component names, click row **5** in the Components table. The information in each field is not saved until you click outside the field.
3. Select **Paroxetine** and modify the levels for this component as shown in the following table.

Cal level	Amount (ng)
Cal_01	25
Cal_02	50
Cal_03	75
Cal_04	100
Cal_05	250
Cal_06	500
Cal_07	750
Cal_08	1000

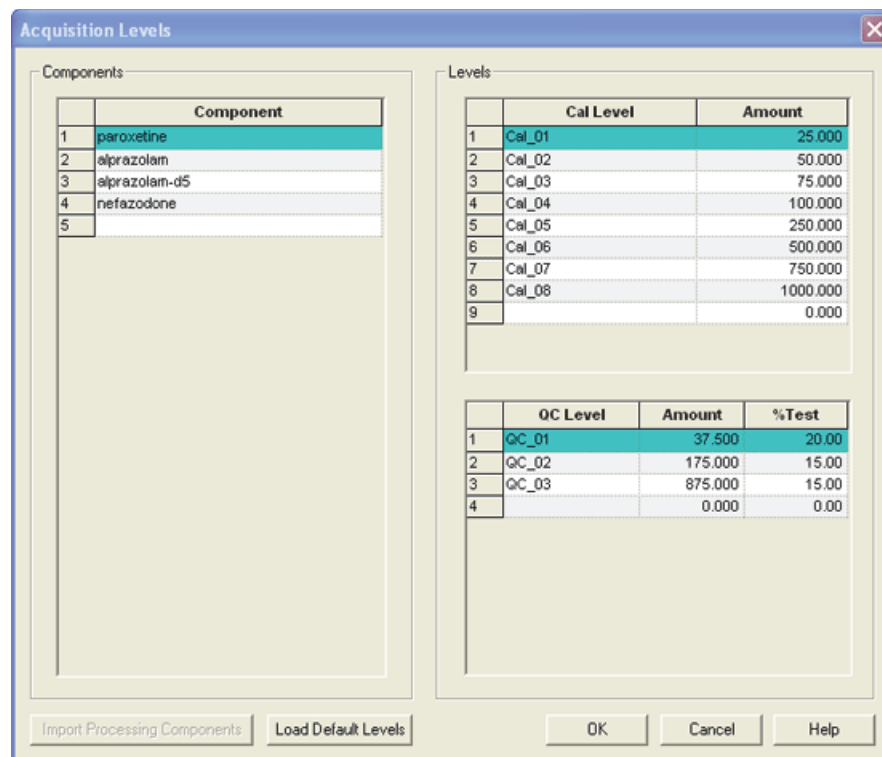
4. To save the new level information for Cal_08, click row **9** in the Levels table.

The information in each field is not saved until you click outside the field.

5. Modify the three QC Level, Amount, and %Test values as shown in the following table.

QC level	Amount	%Test
QC_01	37.5	20
QC_02	175	15
QC_03	875	15

Figure 25. Acquisition Levels dialog box



6. Select **Paroxetine**, right-click anywhere in the Calibration Levels table, and choose **Copy Current Component Levels to All Target Compounds**.

You can see that these same values are now specified for all the components.

7. Click **OK** to save your changes and close the dialog box.
8. Go to [“Exercise 3: Exploring the Data.”](#)

Saving the Workbook and Exiting the LCquan Application

If you want to take a break, you can save your workbook and open it again when you are ready to continue. For instructions about saving a workbook, see [“Saving the Workbook and Exiting the LCquan Application”](#) on [page 59](#).

Exercise 3: Exploring the Data

Use the Explore view to display a multi-component chromatogram for a single sample and experiment with peak detection and integration parameters to see how they affect the chromatogram. In the Explore view, you can determine the optimum parameters used in the processing method. In the Quantitate view, the processing method provides instructions to the LCquan application about how to perform quantitative analysis on raw data.

Opening a Raw Data File and Displaying the Chromatogram

Open a raw data file that represents your data set to determine suitable peak detection and integration parameters. In general, choose a raw data file that corresponds to a low-concentration calibration standard. For this tutorial, you use APN_004.raw.

❖ To enter the Explore view and open the raw data file

1. In the navigation pane, click .

The Create page of the Explore view opens (see [Figure 26](#)). The Chromatogram and Spectrum panes are blank.

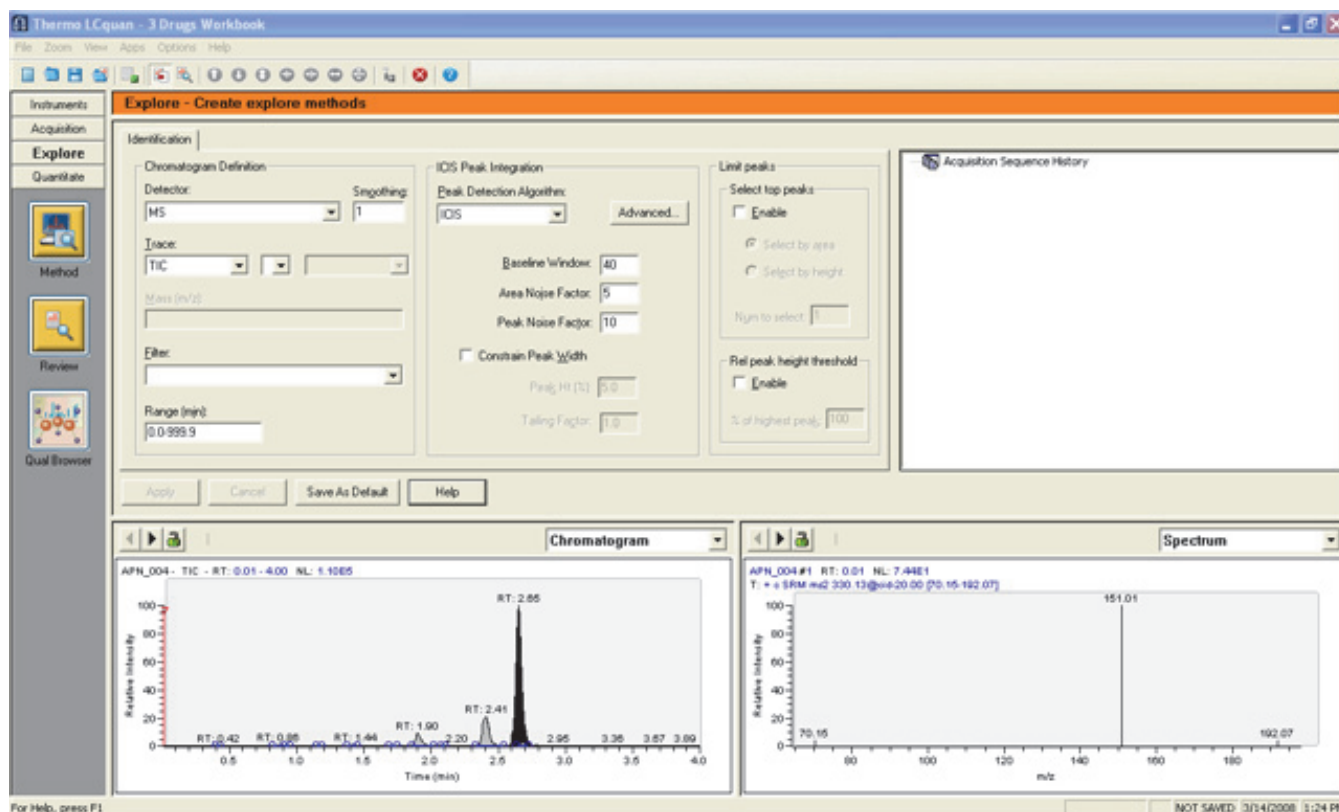
2. Open APN_004.raw:
 - a. From the main menu, choose **File > Open Raw File** to open the Select Rawfile dialog box.
 - b. Browse to the Rawfiles folder:
...\\Xcalibur\\QuanRoot\\Practice Study\\3-Drugs Workbook\\Rawfiles
 - c. Select **APN_004** and click **Open**.

The LCquan application imports the raw data for the APN_004 sample and displays the raw chromatogram on the Create page (see [Figure 26](#)).

Three peaks, which correspond to paroxetine, nefazodone, and both forms of alprazolam, appear in the chromatogram at retention times of 1.90, 2.41, and 2.65 minutes.

The following Create page shows the raw chromatogram for APN_004.

Figure 26. Explore – Create page



Exploring Peak Detection and Integration Parameters

When the LCQuan application acquires data, it creates unique scan filters (see [Table 1](#)) according to the type of experiment you specify in the Instrument Method. When you load a raw data file, the LCQuan application loads the scan filters associated with the raw data file.

Scan filter names are more easily identified when you know the parent ion mass value of the component.

Table 1. Scan filters

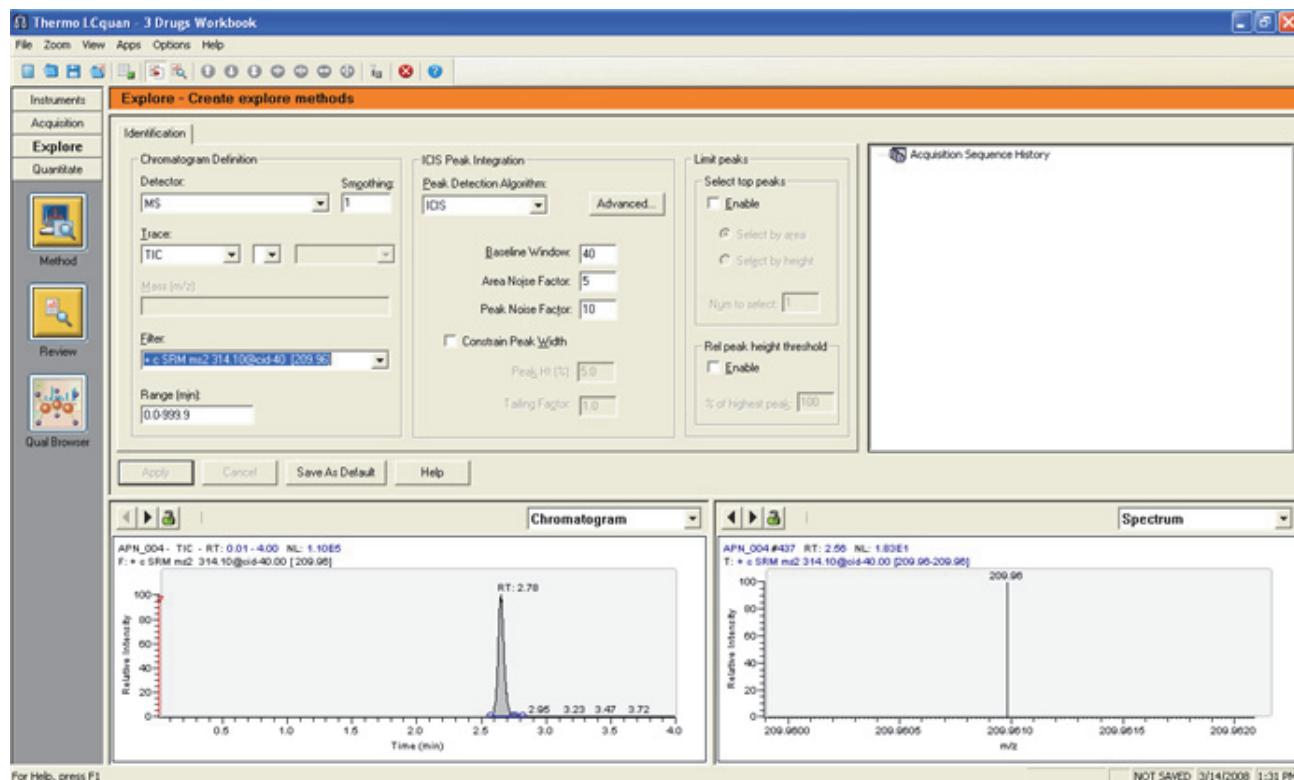
Component name	Parent ion mass <i>m/z</i>	Scan filter
paroxetine	330	+ c SRM ms2 330.13 @cid-20 [70.15-192.07]
nefazodone	470	+ c SRM ms2 470.18 @cid-26 [180.07-274.08]
alprazolam	309	+ c SRM ms2 309.07 @cid-42 [205.02]
alprazolam-d5	314	+ c SRM ms2 314.10 @cid40.00 [209.96]

❖ To relate the chromatogram peaks with the components

1. From the Filter list, select the scan filter for alprazolam_d5 (parent ion mass m/z 314), + c SRM ms2 314.10@cid40.00 [209.96], and click **Apply** (see Figure 27).

On the following Create page, the LCquan application displays the filtered chromatogram for alprazolam-d5. Alprazolam-d5 corresponds to the peak at retention time 2.78 minutes.

Figure 27. Filtered chromatogram for alprazolam-d5 on the Create page



2. Select the scan filter for paroxetine (parent ion mass m/z 330), + c SRM ms2 330.13@cid-20 [70.15-192.07], and click **Apply** to display the filtered chromatogram.

Paroxetine corresponds to the peak at retention time 1.90 minutes.

3. Select the scan filter for nefazodone (parent ion mass m/z 470), + c SRM ms2 470.18@cid-26 [180.07-274.08], and click **Apply** to display the filtered chromatogram.

Nefazodone corresponds to the peak at retention time 2.41 minutes.

4. Select the scan filter for alprazolam (parent ion mass m/z 309), + c SRM ms2 309.07@cid-42 [205.02], and click **Apply** to display the filtered chromatogram.

Alprazolam corresponds to the peak at retention time 2.68 minutes.

5. Select Smoothing: 7 and Peak Detection Algorithm: **Genesis**. Click **Apply**.

Notice the effect on the alprazolam peak.

Note Use the Explore view to experiment with the peak detection and peak integration parameters. Refer to LCquan Help for descriptions of the peak detection and peak integrating parameters.

6. To display the chromatogram for all three components, clear the Filter box and click **Apply**.
7. Click the **Review** icon.

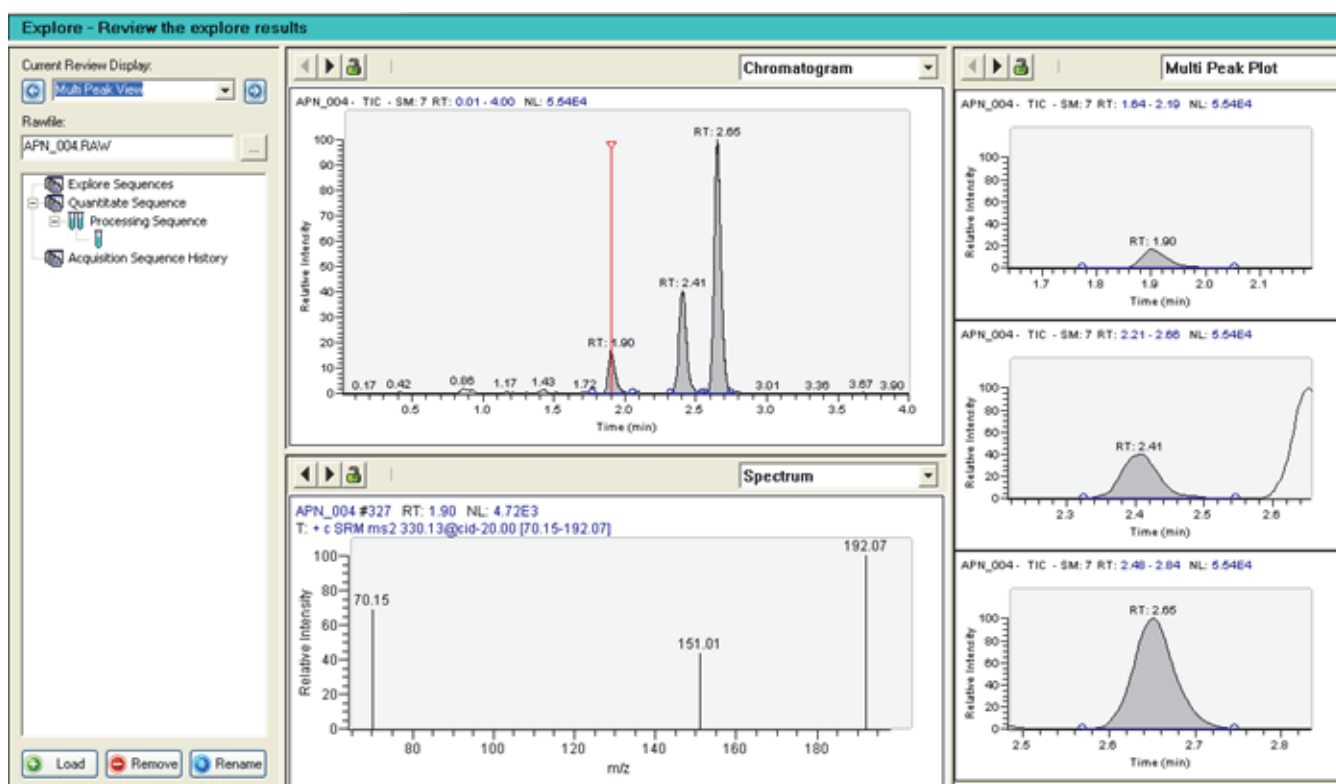
The LCquan application displays the Review page in the Explore view.

Reviewing the Explore Results and Determining Major Product Ions

The Review page displays multiple views of the data. The page shows a chromatogram pane that displays all three peaks, a multi-peak plot pane that individually displays the three peaks, and a spectrum pane that displays the mass-to-charge ratios for the selected peak.

Figure 28 shows the unfiltered chromatogram for the three-drugs data.

Figure 28. Unfiltered chromatogram for the three-drugs data on the Review page



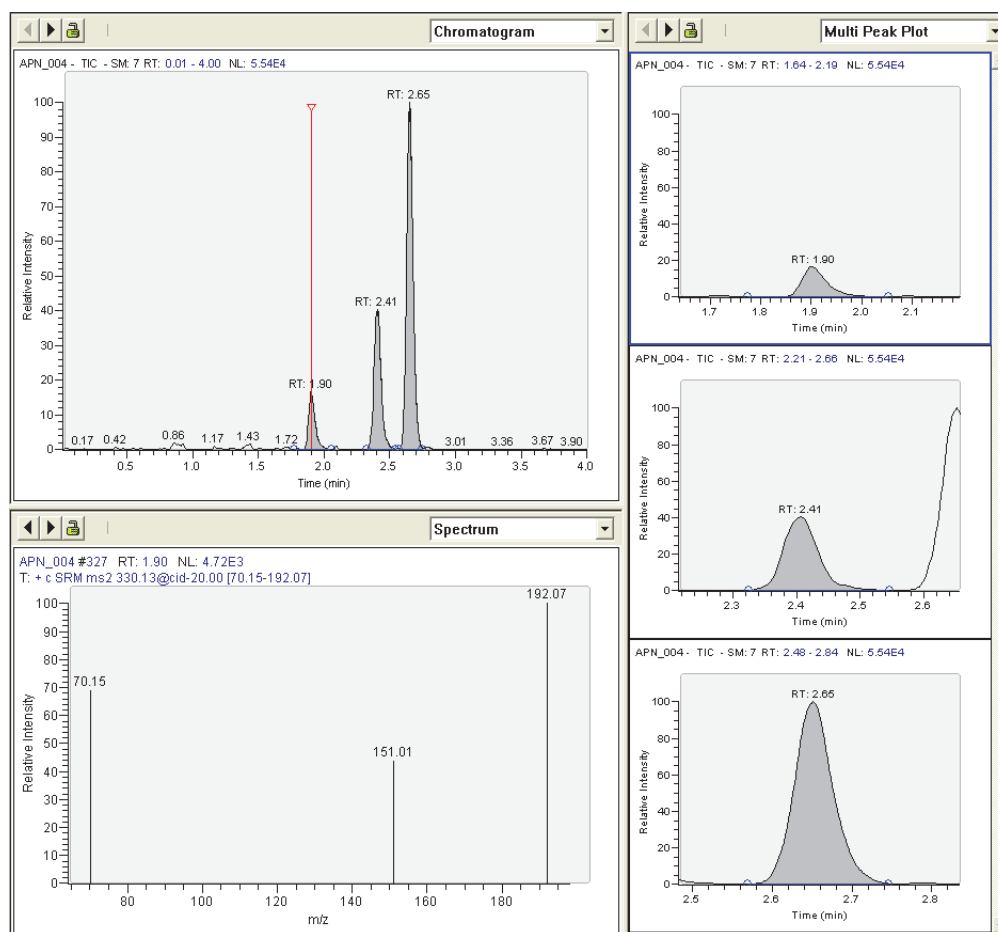
❖ To review the filtered chromatograms

1. In the right-side pane, select **Multi Peak Plot**.
2. Click the top peak (paroxetine) in the Multi Peak Plot.

The Spectrum pane displays the product ion mass spectrum at the maximum intensity.

3. Notice the mass-to-charge ratios (m/z) of the first several ions of highest intensity in the paroxetine mass spectrum (70.15, 151.01, 192.07) (see Figure 29).

Figure 29. Paroxetine data

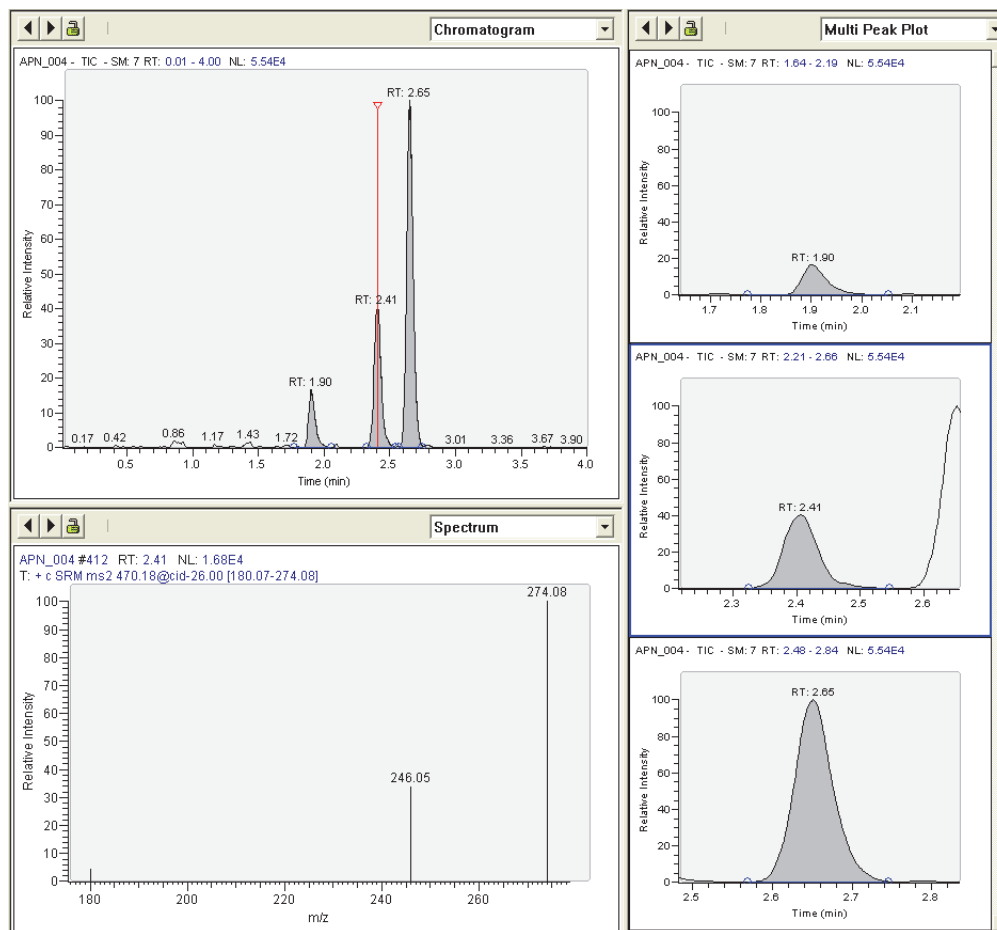


4. Click the second chromatogram peak (nefazodone) in the Multi Peak Plot.

The Spectrum pane displays the product ion mass spectrum at the maximum intensity.

- Notice the mass-to-charge ratios (m/z) of the first several ions of highest intensity in the nefazodone mass spectrum (see [Figure 30](#)).

Figure 30. Nefazodone data



- Click the third chromatogram peak (alprazolam) in the Multi Peak Plot.
The Spectrum pane displays the product ion mass spectrum at the maximum intensity.
- Notice the mass-to-charge ratios (m/z) of the first several ions of highest intensity in the alprazolam mass spectrum (see [Figure 31](#)).

Figure 31. Alprazolam data

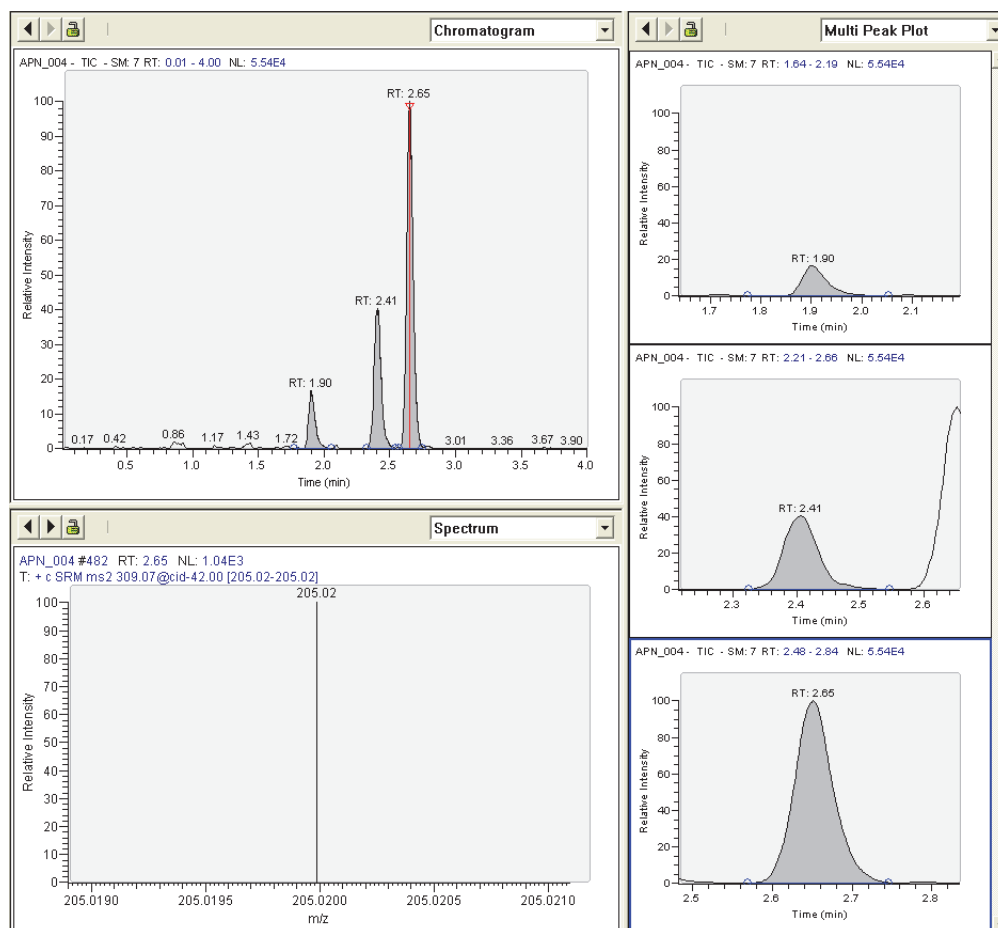


Table 2 lists the mass-to-charge ratios of the three most intense product ions of the three-drugs mass spectra.

Table 2. Mass-to-charge ratios

Paroxetine	Nefazodone	Alprazolam
192.07	274.08	205.02
70.15	246.05	
151.01		

You now have enough information to perform quantitative analysis of the drugs.

- Go to “Exercise 4: Creating a Processing Method.”

Saving the Workbook and Exiting the LCquan Application

If you want to take a break, you can save your workbook and open it again when you are ready to continue. For instructions about saving a workbook, see “Saving the Workbook and Exiting the LCquan Application” on page 59.

Exercise 4: Creating a Processing Method

A processing method provides instructions to the LCQuan application for performing quantitative analysis on raw data. A processing method contains component identification, detection, integration, and calibration information. To create a processing method, use the Create Method pages in the Quantitate view.

Opening the Create Method Page and Using the New Method Wizard

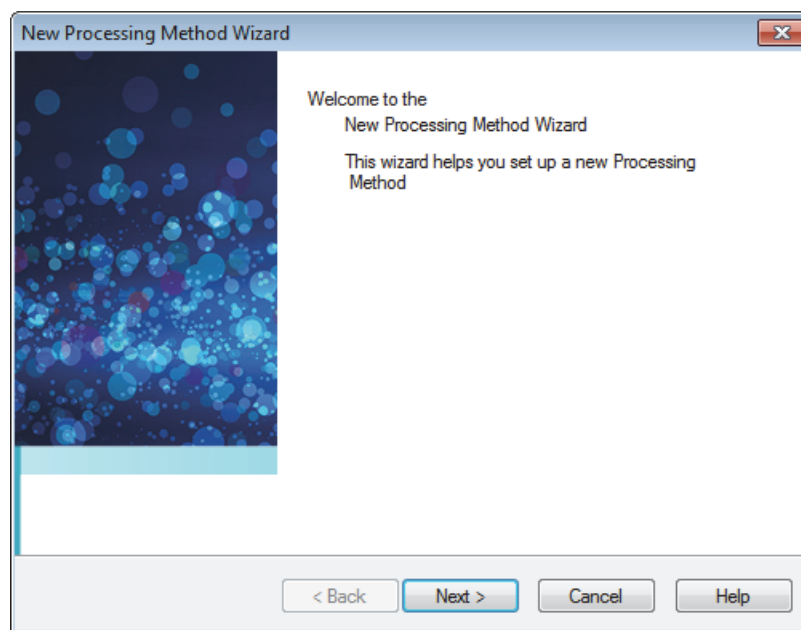
In this exercise, open the Create Method page in the Quantitate view to construct a processing method. When you open this page, the New Processing Method Wizard opens and leads you through the initial steps of creating a method.

❖ **To open the Create Method page of the Quantitate view and start the New Processing Method Wizard**

1. From any view, click  in the LCQuan navigation pane.

The New Processing Method Wizard opens (see [Figure 32](#)).

Figure 32. Welcome page of the New Method Wizard



2. Click **Next**, and then select the **Create New Method** option and the **Initialize with Acquisition Component Names** check box (see [Figure 33](#)).

Figure 33. Create new method options page

Create a new empty method or import a method from legacy data files.

☒ Create new method
☒ Initialize with acquisition component names
☐ Import existing method

3. Click **Next**, and then select the **Calibrate Using Internal Standards** option to specify that you calibrate with an internal standard (see [Figure 34](#)).

Figure 34. Calibrate options page

Select the type of standards to use for the new method.

☒ Calibrate using Internal Standards
☐ Calibrate using External Standards

4. Click **Next**.
5. Open a raw data file from your data set to determine the peak detection and integration parameters, as follows:
 - a. Click **Browse** to find the APN_004.raw data file in the Rawfiles folder in your new workbook.

The example uses this default path:

...\\Xcalibur\\QuanRoot\\Practice Study\\3-Drugs Workbook\\Rawfiles\\APN_04

- b. Click **Open** and click **Next** (see [Figure 35](#)).

Figure 35. Raw data files page

Example raw file used during method development to view the effects caused by changes in detection and integration parameters

Raw File

Example Study\\3-Drugs Workbook\\Rawfiles\\APN_015.RAW

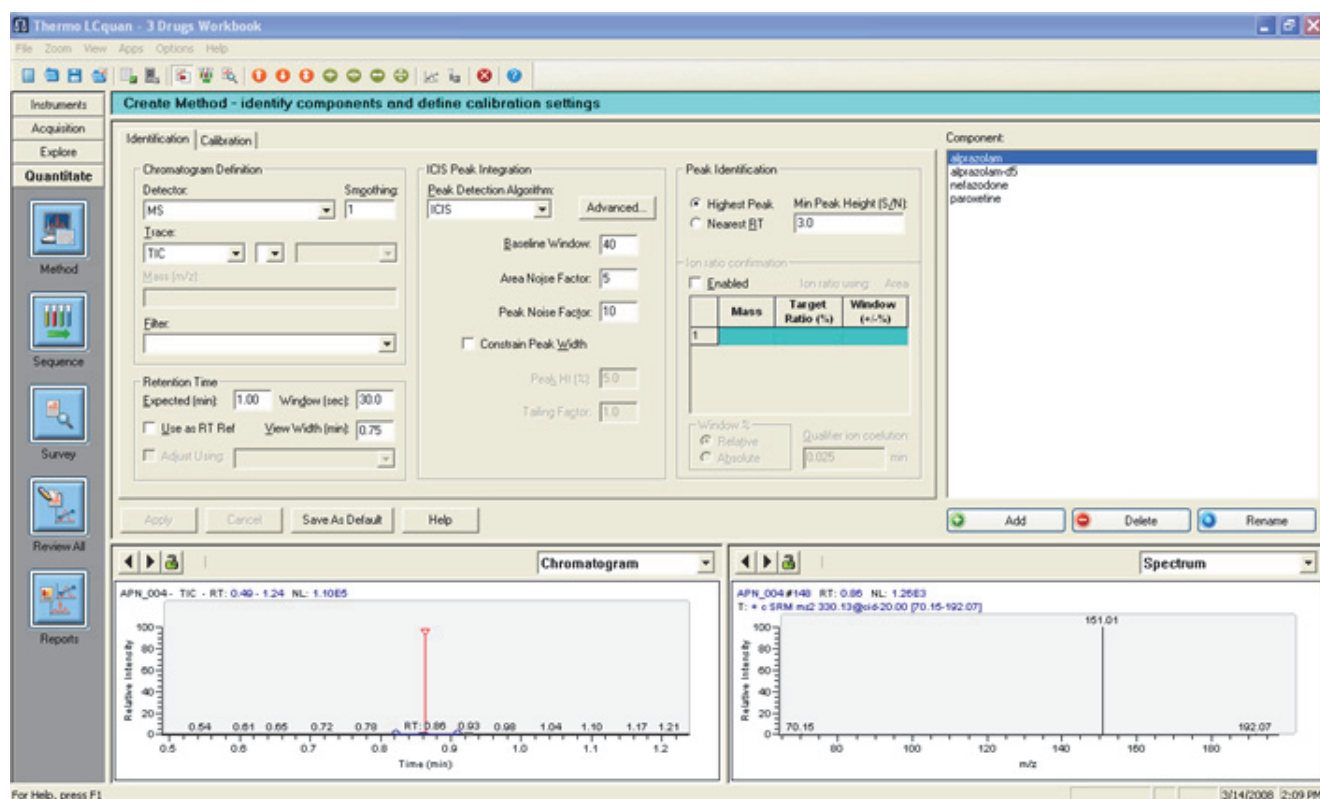
Browse...

Tip In general, choose a raw data file that corresponds to a low-concentration calibration standard.

- Click **Finish** to exit the New Method Wizard.

The Create Method page opens (see Figure 36) for you to begin building the processing method.

Figure 36. Create Method page in the Quantitate view



Identifying the Internal Standard

In the last exercise, you imported the component names and calibration levels from the acquisition sequence into the processing method. In this exercise, you will identify the internal standard.

❖ To identify alprazolam-d5 as the internal standard compound and specify its calibration parameters

- On the Create Method page in the Quantitate view, select **Alprazolam-d5** in the Component list.
- Click the **Calibration** tab to display the calibration settings.
- Enter the values and options as indicated in Figure 37.

Figure 37. Calibration page showing the selection of Alprazolam-d5

Identification Calibration

Component Type

☐ Target Compound

☒ ISTD

Target Units:

ISTD Amount:

1.000

ISTD Units:

ng/ml

Target Compound

ISTD:

alprazolam-d5

Isotope %...

Calibration Curve Type:

Average RF

Response

☒ Area

☐ Height

Weighting

☒ Equal

☐ 1/X

☐ 1/X^2

☐ 1/Y

☐ 1/Y^2

☐ 1/s^2

Origin

☒ Ignore

☐ Force

☐ Include

4. Click **Apply** to accept the settings.

If necessary, click **OK** to dismiss the warning box.

Specifying Calibration Curve Parameters for the Target Compounds

❖ To specify the calibration curve parameters for the target compounds

1. On the Create Method page in the Quantitate view, specify the calibration curve parameters for paroxetine:
 - a. In the Component list, select **Paroxetine**.
 - b. On the Calibration page, enter the values and options as indicated in [Figure 38](#).

Figure 38. Calibration page

Identification Calibration

Component Type

☒ Target Compound

☐ ISTD

Target Units:

ng/ml

ISTD Amount:

1.000

ISTD Units:

Target Compound

ISTD:

alprazolam-d5

Isotope %...

Calibration Curve Type:

Linear

Response

☒ Area

☐ Height

Weighting

☒ Equal

☐ 1/X

☐ 1/X^2

☐ 1/Y

☐ 1/Y^2

☐ 1/s^2

Origin

☐ Ignore

☐ Force

☒ Include

- c. Click **Apply** to apply the settings.
2. On the Create Method page in the Quantitate view, specify the calibration curve parameters for nefazodone, as follows:
 - a. In the Component list, select **Nefazodone**.
 - b. On the Calibration page, enter the values and options as indicated in [Figure 39](#).

Figure 39. Calibration page

The screenshot shows the 'Calibration' tab of a software interface. The 'Component Type' section on the left has 'Target Compound' selected. The 'Target Units' field contains 'ng/ml'. The 'ISTD Amount' field contains '1.000'. The 'ISTD Units' field is empty. The 'Target Compound' field on the right contains 'alprazolam-d5'. The 'Isotope %...' field is empty. The 'Calibration Curve Type' dropdown is set to 'Linear'. The 'Response' section has 'Area' selected. The 'Weighting' section has '1/X' selected. The 'Origin' section has 'Ignore' selected.

- c. Click **Apply** to apply the settings.
3. On the Create Method page in the Quantitate view, specify the calibration curve parameters for alprazolam:
 - a. Select **Alprazolam** in the Component box.
 - b. On the Calibration page, enter the values and options indicated in [Figure 40](#).

1 Tutorial Exercises

Exercise 4: Creating a Processing Method

Figure 40. Calibration page

The screenshot shows the 'Calibration' tab of a software interface. It contains several configuration options for a target compound. On the left, under 'Component Type', 'Target Compound' is selected. Below this, 'Target Units' is set to 'pg/ml', 'ISTD Amount' is '1.000', and 'ISTD Units' is empty. On the right, 'Target Compound' is 'alprazolam-d5'. Below this is an 'Isotope %...' button. 'Calibration Curve Type' is set to 'Linear'. Under 'Response', 'Area' is selected. Under 'Weighting', '1/X' is selected. Under 'Origin', 'Ignore' is selected.

- c. Click **Apply** to apply the settings.
4. Go to the next topic: [“Entering Component Identification Parameters and Integrating the Peaks.”](#)

Entering Component Identification Parameters and Integrating the Peaks

When the LCquan application acquires data, it creates unique scan filters according to the type of experiment you specify in the Instrument Method. To quantitate paroxetine, nefazodone, and alprazolam, you must filter the chromatogram.

❖ **To specify the identification settings that you determined in the Explore view for each of the drugs**

1. Click the **Identification** tab (see [Figure 41](#)).

Figure 41. Identification page

The screenshot shows the 'Identification' tab of the LCQuan software interface. It is divided into two main sections: 'Chromatogram Definition' and 'ICIS Peak Integration'.

Chromatogram Definition:

- Detector:** MS
- Smoothing:** 1
- Trace:** Mass Range
- Mass (m/z):** 209.96
- Filter:** + c SRM ms2 314.10@cid-40.00 [209.96-209.9]
- Retention Time:**
 - Expected (min):** 1.00
 - Window (sec):** 30.0
 - ☒ **Use as RT Ref**
 - View Width (min):** 0.75
 - ☐ **Adjust Using:** (empty dropdown)

ICIS Peak Integration:

- Peak Detection Algorithm:** ICIS
- Baseline Window:** 30
- Area Noise Factor:** 5
- Peak Noise Factor:** 10
- ☐ **Constrain Peak Width**
- Peak Ht (%):** 5.0
- Tailing Factor:** 1.0

2. In the Component list, select **Alprazolam-d5**.
3. Enter the settings for Alprazolam-d5, as follows:
 - In the Chromatogram Definition area:
 - Detector: **MS**
 - Smoothing: **1**
 - Trace: **Mass Range** (specifies a mass range, not TIC, chromatogram)
 - Mass: **209.96** (important mass that you determined in the Explore view)
 - Filter: **+ c SRM ms2 314.10@cid-40 [209.98]** (scan filter for m/z 314 parent mass)
 - In the ICIS Peak Integration area:
 - Peak Detection Algorithm: **ICIS**
 - Baseline Window: **30**
 - Area Noise Factor: **5**
 - Peak Noise Factor: **10**

Click **Apply** to apply the settings.

The LCQuan application applies the settings to the data in the raw data file and automatically displays the resulting filtered chromatogram in the Chromatogram pane.

4. Specify the expected retention time of alprazolam-d5, as follows (see [Figure 42](#)):

Figure 42. Identification page showing the component retention time for Alprazolam-d5

The screenshot shows the 'Identification' tab of the LCQuan software interface. It is divided into two main sections: 'Chromatogram Definition' and 'Retention Time'.
 In the 'Chromatogram Definition' section:
 - 'Detector' is set to 'MS'.
 - 'Smoothing' is set to '1'.
 - 'Trace' is set to 'Mass Range'.
 - 'Mass (m/z)' is set to '209.96'.
 - 'Filter' is set to '+ c SRM ms2 314.10@cid-40.00 [209.96]'.
 In the 'Retention Time' section:
 - 'Expected (min)' is set to '2.65'.
 - 'Window (sec)' is set to '30.0'.
 - The 'Use as RT Ref' checkbox is checked.
 - 'View Width (min)' is set to '0.75'.
 - The 'Adjust Using' dropdown is empty.

- a. Inspect the mass chromatogram view and notice the retention time of alprazolam-d5.
 Each component peak appears with the appropriate retention time label. In this example the alprazolam-d5 peak is labeled RT:2.65.
- b. In the Expected (min) text box, type **2.65** to specify the retention time of alprazolam-d5 as 2.65 minutes.
- c. In the Window box, type **30.0** to specify the peak detection time search window as 30 seconds wide (centered about the retention time).
- d. Select the **Use as RT Ref** check box.
- e. In the View Width box, type **0.75** to specify a chromatogram view width of 0.75 minutes.
- f. Click **Apply** to save the component identification information for alprazolam-d5.

The LCQuan application automatically integrates the peak according to the specified parameters. The integrated portion of the peak appears gray, and blue identification markers appear at the starting and ending points of the peak integration. The blue line that connects the integration markers indicates the baseline. See [Create Method page on page 37](#).

5. In the Component list, select **Paroxetine**.

6. Enter the settings for paroxetine, as follows (see [Figure 43](#)):

Figure 43. Identification page for paroxetine

The screenshot shows the 'Identification' tab of a software interface. It is divided into two main panels. The left panel, 'Chromatogram Definition', contains settings for the detector (MS), smoothing (1), trace (Mass Range), mass (m/z) (192.07), and filter (+ c SRM ms2 330.13@cid-20.00 [70.15-192.0]). The right panel, 'ICIS Peak Integration', contains settings for the peak detection algorithm (ICIS), baseline window (20), area noise factor (5), peak noise factor (10), and options for constraining peak width, peak height (5.0%), and tailing factor (1.0). There are also retention time settings at the bottom left.

- In the Chromatogram Definition area:
 - Detector: **MS**
 - Smoothing: **1**
 - Trace: **Mass Range**
 - Mass (*m/z*): **192.07**
 - Filter: **+ c SRM ms2 330.13@cid-20 [70.15-192.07]**
- In the ICIS Peak Integration area:
 - Peak Detection Algorithm: **ICIS**
 - Baseline Window: **20**
 - Area Noise Factor: **5**
 - Peak Noise Factor: **10**

Click **Apply** to apply the settings.

1 Tutorial Exercises

Exercise 4: Creating a Processing Method

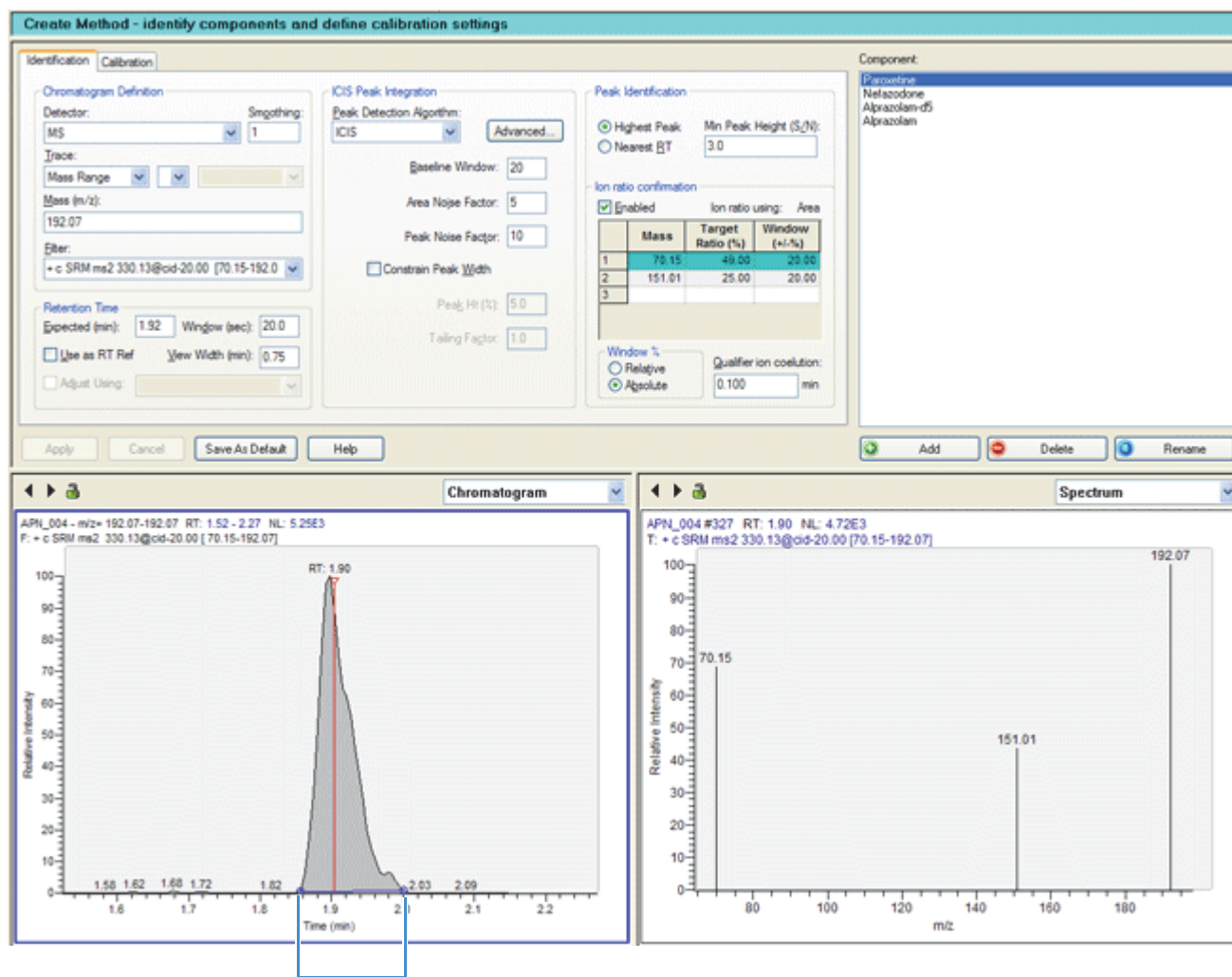
7. Specify the expected retention time of paroxetine, as follows (see Figure 44):

Figure 44. Identification page showing the component retention time for paroxetine

- Inspect the mass chromatogram and note the retention time of paroxetine. In this example the paroxetine peak is labeled *RT:1.9*.
 - In the Expected (min) text box, type **1.9** to specify the retention time of paroxetine as 1.9 minutes.
 - In the Ion ratio confirmation area, select the **Enabled** check box.
 - Specify the following settings for the mass, target ratio, and window options.
- | Mass | Target ratio | Window |
|--------|--------------|--------|
| 70.15 | 49 | 20 |
| 151.01 | 25 | 20 |
- Select the **Absolute** option for the Window% setting.
 - Set the Qualifier ion coelution to **0.1** minutes.
8. Integrate the paroxetine chromatogram peak:
- Click the chromatogram peak in the Chromatogram pane.
 - Select the **Adjust Using: Alprazolam-d5** check box to use the retention time of alprazolam-d5 to adjust for the retention time drift of paroxetine.
 - Click **Apply** to integrate the chromatogram peak.

Figure 45 shows the integrated mass chromatogram for paroxetine.

Figure 45. Integrated mass chromatogram for paroxetine on the Create Method page



9. In the Component list, select **Nefazodone**.

10. Specify the settings for nefazodone, as follows (see Figure 46):

Figure 46. Identification page for nefazodone

The screenshot shows the 'Identification' tab of the LCQuan software interface. The 'Chromatogram Definition' section is configured with the following settings: Detector: MS, Smoothing: 1, Trace: Mass Range, Mass (m/z): 274.08, and Filter: + c SRM ms2 470.18@cid-26.00 [180.07-274.08]. The 'ICIS Peak Integration' section is configured with: Peak Detection Algorithm: ICIS, Baseline Window: 20, Area Noise Factor: 5, Peak Noise Factor: 10, Constrain Peak Width: unchecked, Peak Ht (%): 5.0, and Tailing Factor: 1.0. The 'Retention Time' section includes Expected (min): 1.00, Window (sec): 30.0, Use as RT Ref: unchecked, View Width (min): 0.75, and Adjust Using: empty.

- In the Chromatogram Definition area:
 - Detector: **MS**
 - Smoothing: **1**
 - Trace: **Mass Range**
 - Mass (*m/z*): **274.08**
 - Filter: **+ c SRM ms2 470.18@cid-26.00 [180.07-274.08]**
- In the ICIS Peak Integration area:
 - Peak Detection Algorithm: **ICIS**
 - Baseline Window: **20**
 - Area Noise Factor: **5**
 - Peak Noise Factor: **10**

Click **Apply** to apply the settings.

11. Specify the expected retention time of nefazodone, as follows (see Figure 47):

Figure 47. Identification page showing the component retention time for nefazodone

Identification | Calibration

Chromatogram Definition

Detector: MS Smoothing: 1

Trace: Mass Range

Mass (m/z): 274.08

Filter: + c SRM ms2 470.18@cid-26.00 [180.07-274.]

Retention Time

Expected (min): 2.41 Window (sec): 30.0

☐ Use as RT Ref View Width (min): 0.75

☐ Adjust Using:

ICIS Peak Integration

Peak Detection Algorithm: ICIS Advanced...

Baseline Window: 20

Area Noise Factor: 5

Peak Noise Factor: 10

☐ Constrain Peak Width

Peak Ht (%): 5.0

Tailing Factor: 1.0

Peak Identification

☒ Highest Peak Min Peak Height (S/N): 3.0

☐ Nearest RT

☒ Ion ratio confirmation Enabled Ion ratio using: Area

	Mass	Target Ratio (%)	Window (+/-%)
1	246.05	30.00	20.00
2	180.07	8.00	20.00
3			

Window % ☐ Relative ☒ Absolute

Qualifier ion coelution: 0.100 min

- Inspect the mass chromatogram and notice the retention time of paroxetine.

In this example, the nefazodone peak is labeled 2.41.

- In the Expected (min) text box, type **2.41** to specify the retention time of nefazodone as 2.41 min.
- In the Ion ratio confirmation area, select the **Enabled** check box.
- Enter these settings for the mass, target ratio, and window options.

Mass	Target ratio	Window
246.05	30	20
180.07	8	20

- Select the **Absolute** option for the Window% setting.
- Set the Qualifier ion coelution to **0.1** minutes.

12. Integrate the nefazodone chromatogram peak, as follows:

- Click the chromatogram peak (see Figure 48) in the Chromatogram pane.
- Click **Apply** to integrate the chromatogram peak (see Figure 49).

1 Tutorial Exercises

Exercise 4: Creating a Processing Method

Figure 48. IRC chromatogram peaks for nefazodone

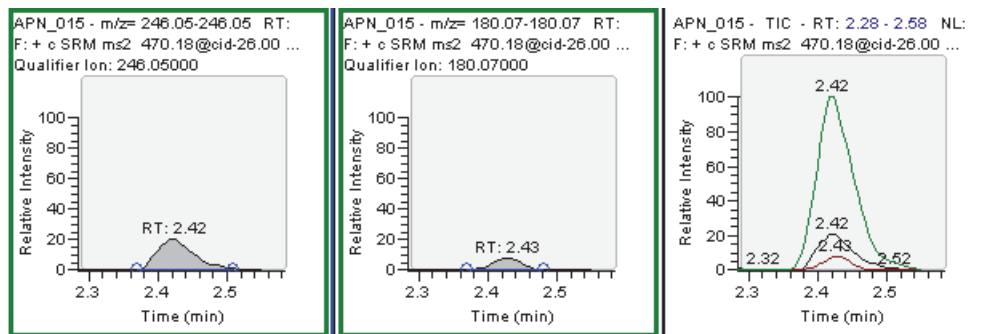
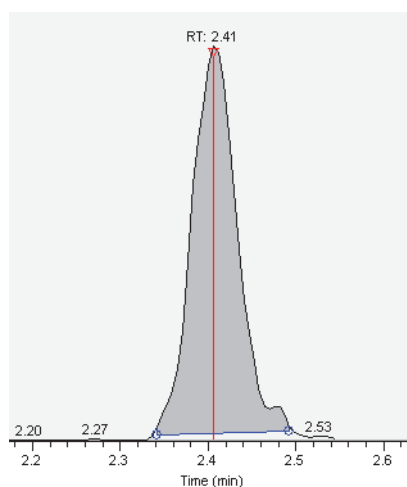


Figure 49. Integrated nefazodone peak



13. In the Component list, select **Alprazolam** and specify these settings on the Identification page in the Chromatogram Definition area (see [Figure 50](#)):

- Detector: **MS**
- Smoothing: **1**
- Trace: **Mass Range**
- Mass (m/z): **205.02**
- Filter: **+ c SRM ms2 309.07@cid-42.00 [205.02]**
- In the ICIS Peak Integration area:
- Peak Detection Algorithm: **ICIS**
- Baseline Window: **20**
- Area Noise Factor: **5**
- Peak Noise Factor: **10**

Figure 50. Identification page for alprazolam

The screenshot shows the 'Identification' tab of a software interface. It is divided into two main sections: 'Chromatogram Definition' and 'ICIS Peak Integration'.

Chromatogram Definition:

- Detector:** MS
- Smoothing:** 1
- Trace:** Mass Range
- Mass (m/z):** 205.02
- Filter:** + c SRM ms2 309.07@cid-42.00 [205.02]
- Retention Time:**
 - Expected (min):** 1.00
 - Window (sec):** 30.0
 - ☐ Use as RT Ref
 - View Width (min):** 0.75
 - ☐ Adjust Using:

ICIS Peak Integration:

- Peak Detection Algorithm:** ICIS
- Baseline Window:** 20
- Area Noise Factor:** 5
- Peak Noise Factor:** 10
- ☐ Constrain Peak Width
- Peak Ht (%):** 5.0
- Tailing Factor:** 1.0

14. Click **Apply** to apply the settings.
15. Specify the expected retention time of alprazolam:
 - a. Inspect the mass chromatogram and notice the retention time of alprazolam.
In the following example, the alprazolam peak is labeled 2.68.
 - b. In the Expected (min) text box, type **2.68** to specify the retention time of alprazolam as 2.68 min.
16. Integrate the alprazolam chromatogram peak, as follows:
 - a. Click the chromatogram peak in the Chromatogram pane.
 - b. Click **Apply** to integrate the chromatogram peak (Figure 51).

Figure 51. Integrated alprazolam peak



Saving the Processing Method

When you have finished creating the processing method, choose **File > Save** to save the settings in the workbook and go to [“Exercise 5: Creating a Processing Sequence.”](#)

Saving the Workbook and Exiting the LCquan Application

If you want to take a break, you can save your workbook and open it again when you are ready to continue. For instructions about saving a workbook, see [“Saving the Workbook and Exiting the LCquan Application”](#) on [page 59](#).

Exercise 5: Creating a Processing Sequence

This exercise creates a processing sequence by using the Create Sequence pages of the Quantitate view. A processing sequence provides instructions to the LCquan application about how to process a batch of raw data files. It consists of a list of sample data files and includes information about sample type and calibration or quality control levels.

❖ To create a new processing sequence for the three-drugs example

1. Open a Create Sequence page and use the New Processing Sequence Wizard.
2. Associate raw data files with sample types.
3. Review the processing sequence.

Opening the Create Sequence Page and Using the New Processing Sequence Wizard

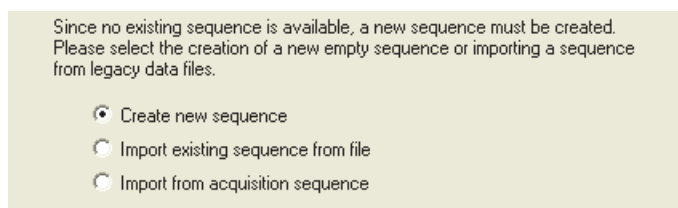
In this exercise, use the Create Sequence pages to create a new processing sequence. When you click the Sequence icon, the LCquan application opens the New Processing Method Wizard unless you have not previously created a processing sequence. If you have not previously created a processing sequence, LCquan opens the Create Sequence page.

❖ To create a new processing sequence (or import the acquisition sequence) for the three-drugs example

1. Click  and click the **Sequence** icon to open the Welcome page of the New Processing Sequence Wizard.
2. Click **Next**, select the **Create New Sequence** option, and click **Next** again ([Figure 52](#)).

Note To make the processing sequence the same as the acquisition sequence that you created earlier, select Import from Acquisition Sequence, click Next and Finish, and go to [“Processing the Raw Data Files”](#) on [page 50](#).

Figure 52. Create new sequence options page



3. Click **Finish** to exit the New Processing Method Wizard.

The Create Sequence (edit) page opens. You can now begin defining the processing sequence.

Associating Raw Data Files with Sample Types

When you use the LCQuan application to perform quantitative analysis, you usually copy the acquisition sequence from the current workbook to use as the processing sequence, or you import an existing sequence from another workbook or legacy file.

Because this is a tutorial, you do not need to create an acquisition sequence because the raw data already exists. For practice, you created an acquisition sequence in “[Exercise 2: Creating an Acquisition Sequence](#)” on [page 7](#). This acquisition sequence associated raw data files with calibration levels or unknowns, specified that there are four components, and specified the calibration level amounts.

Use the Create Sequence pages to build a processing sequence by associating each raw data file in the data set with one of the following sample types: standard, QC, blank, or unknown; then associate the standards and QCs with the appropriate calibration level or quality control level.

❖ To build a new processing sequence for the three-drugs example

1. Locate the three-drugs raw data set, as follows:
 - a. In the directory tree in the Available Files pane, browse to find the Rawfiles folder in your current workbook. This example uses the following default path:
 ...\\Xcalibur\\QuanRoot\\Practice Study\\3-Drugs Workbook\\Rawfiles
 - b. Click the **Rawfiles** folder.

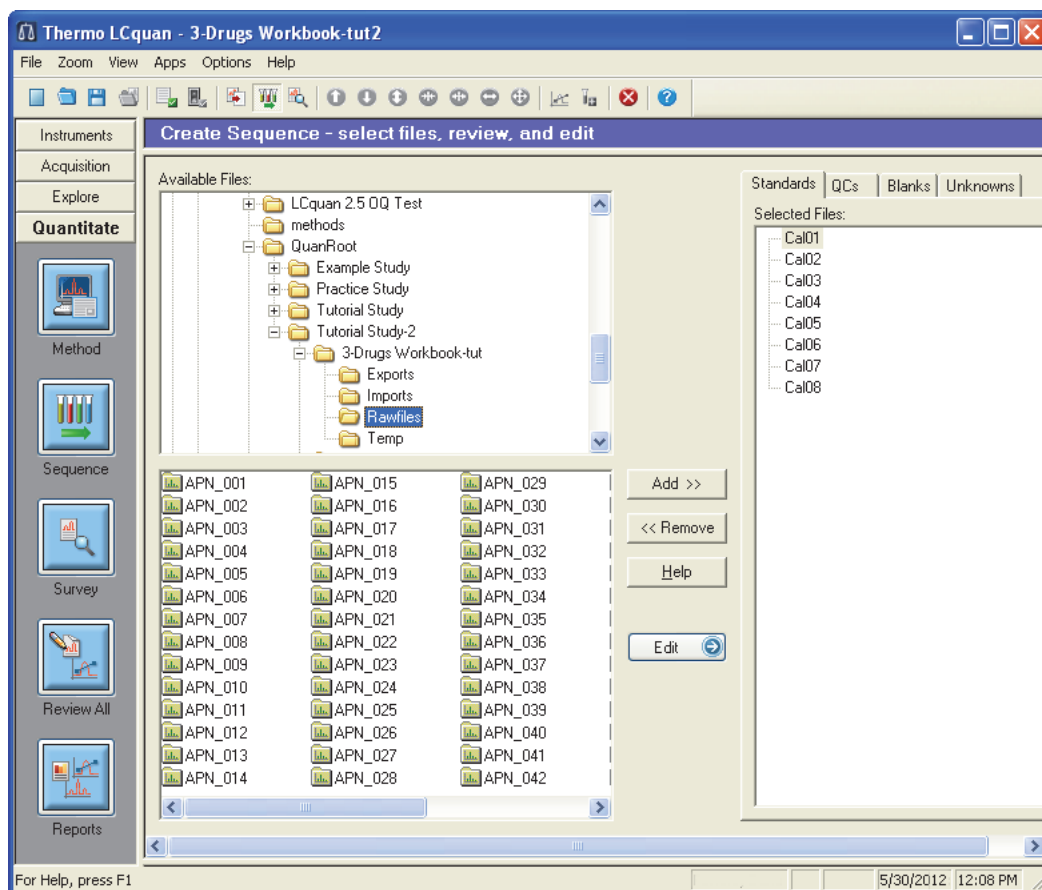
The LCQuan application displays all of the three-drugs raw data files in the Available Files pane.

1 Tutorial Exercises

Exercise 5: Creating a Processing Sequence

Figure 53 shows the raw data files in the three-drugs data set.

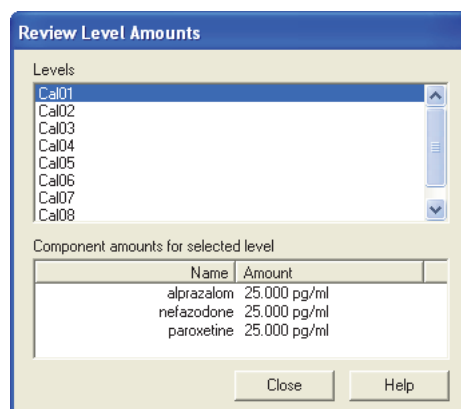
Figure 53. Create Sequence (create) page



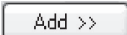
2. Associate the calibration data files with the calibration levels, as follows:
 - a. Click the **Standards** tab to display the calibration levels (Cal01, Cal02, and so on).
 - b. To display the amount of standard sample associated with a calibration level, right-click the calibration level in the Selected Files pane and choose **Show Level Amounts**.

The Review Level Amounts dialog box opens (see [Figure 54](#)).

Figure 54. Review Level Amounts dialog box showing the component amounts



The eight calibration levels contain the amounts of each drug that you specified when you set the acquisition levels.

- c. Click **Close** in the Review Level Amounts dialog box.
- d. In the Selected Files pane, click **Cal01** to select the first calibration level.
- e. In the Available Files pane, click **APN_003** to select the file.
- f. Click .

The LCQuan application adds the selected raw data file to the Cal01 calibration level in the Selected Files pane.

Tip You can also add a raw data file to a calibration level file by dragging the raw data file to the appropriate level in the Selected Files pane.

- g. Add the raw data files to the calibration levels as shown in [Table 3](#).

Table 3. Calibration levels

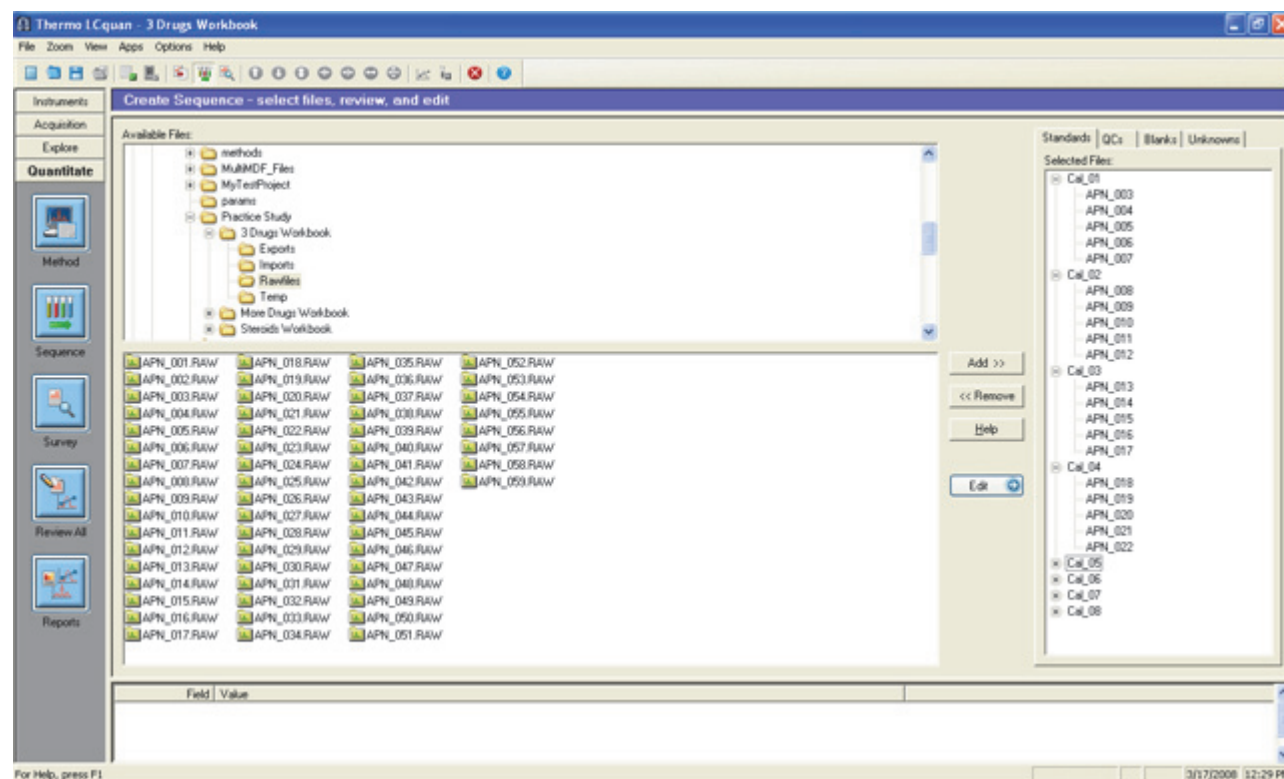
Calibration level	Raw data file
Cal01	APN_003 – APN_007
Cal02	APN_008 – APN_012
Cal03	APN_013 – APN_017
Cal04	APN_018 – APN_022
Cal05	APN_023 – APN_027
Cal06	APN_028 – APN_032
Cal07	APN_033 – APN_037
Cal08	APN_038 – APN_042

1 Tutorial Exercises

Exercise 5: Creating a Processing Sequence

Figure 55 shows the raw data files associated with the calibration levels.

Figure 55. Create Sequence (create) page



Tip If necessary, you can use the sample type and concentration information in the File Information pane at the bottom of the view to identify the content of the datafiles.

3. Associate the data files with the QC levels, as follows:
 - a. Click the **QC** tab in the Selected Files pane.
 - b. In the QCs pane, select the **QC_01** level.
 - c. In the Available Files pane, select the **APN_045** file.
 - d. Click .

The LCQuan application adds the selected raw data file to the QC_01 level in the Selected Files pane.

- e. Associate the raw data files and QC levels as shown in this table.

QCs	Raw data file
QC_01	APN_045 – APN_047
QC_02	APN_048 – APN_050
QC_03	APN_051 – APN_053

4. Associate the data files with the blank sample type, as follows.
 - a. Click the **Blanks** tab in the Selected Files pane.
 - b. In the Available Files pane, hold down SHIFT and select **APN_001** and **APN_002**.
 - c. Click .

The LCQuan application adds the selected raw data files to the Blank pane.

- d. In the Available Files pane, hold down SHIFT and select **APN_043** and **APN_044**.
 - e. Click .

The LCQuan application adds the selected raw data files to the Blank pane.

- f. In the Available Files pane, hold down SHIFT and select **APN_054** and **APN_055**.
 - g. Click .

The LCQuan application adds the selected raw data files to the Blank pane.

5. Associate the remaining raw data files with the unknown sample type, as follows:
 - a. Click the **Unknowns** tab in the Selected Files pane.
 - b. In the Available Files pane, select **APN_056**.
 - c. Hold down SHIFT and select **APN_059** to select all unknowns.
 - d. Click **Add**.

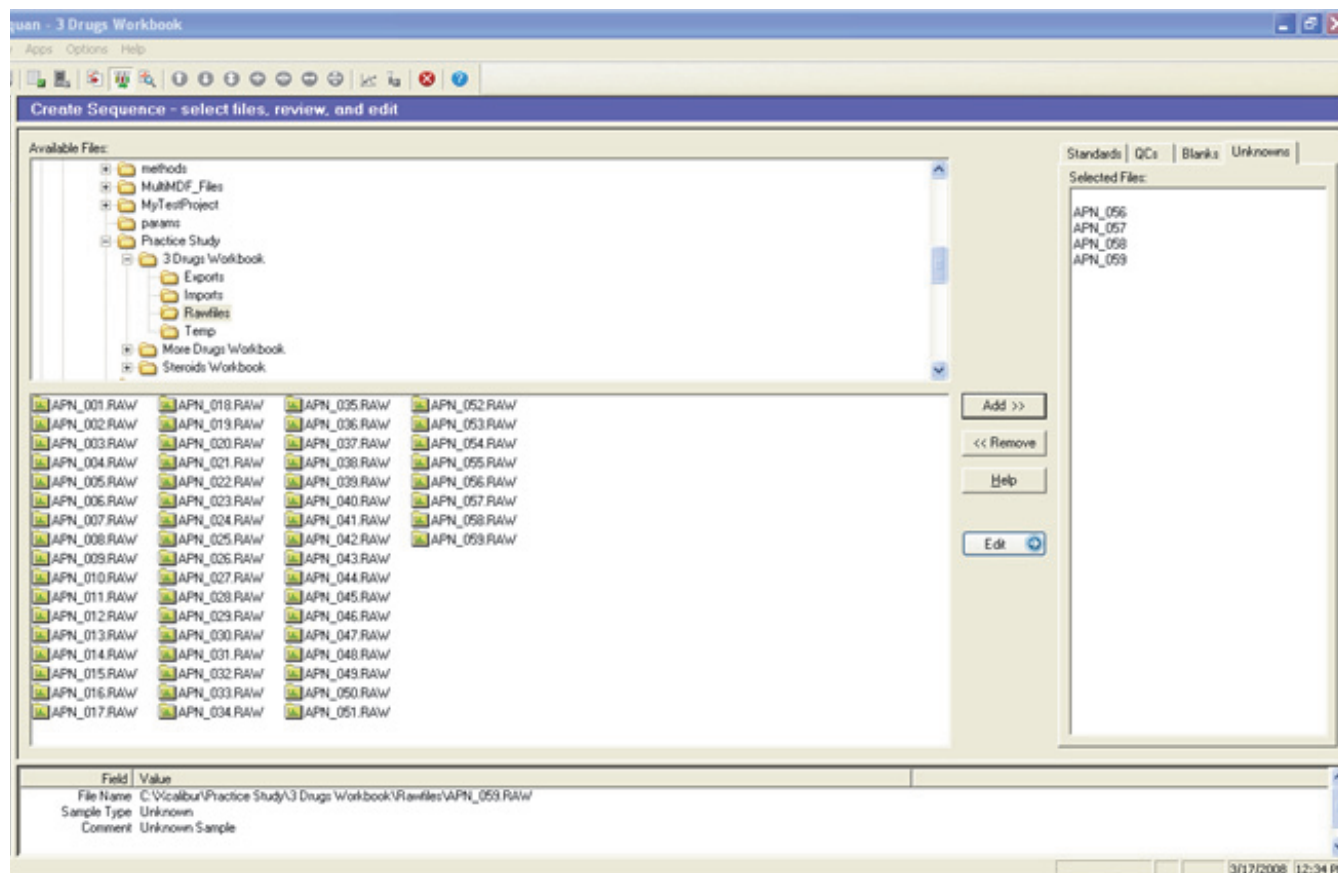
The LCQuan application adds the four selected raw data files to the Unknowns pane.

1 Tutorial Exercises

Exercise 5: Creating a Processing Sequence

Figure 56 shows the proper association of raw data files to the unknown sample type for the three-drugs example.

Figure 56. Create Sequence (create) page



Reviewing the Processing Sequence

After you associate all of the raw data files with sample types, review the processing sequence to ensure it is correct.

❖ To review the processing sequence

1. Click  to open the Create Sequence (edit) page and display the new processing sequence (see Figure 57).

Figure 57. Raw data files for the three-drugs example and their calibration levels

	Sample Type	Filename	Path	Sample ID	Level	ISTD Corr Amt	Dil Factor	Vial Pos	Inj Vol	Sample Vol	Sample Wt
1	Standard	APN_001	C:\Xcalibur\Q	APN002	Cal_01	0.000	1.000	Tray01:2	30.00	0.000	0.000
2	Standard	APN_002	\Rawfiles	APN003	Cal_01	0.000	1.000	Tray01:2	30.00	0.000	0.000
3	Standard	APN_003	\Rawfiles	APN004	Cal_01	0.000	1.000	Tray01:2	30.00	0.000	0.000
4	Standard	APN_004	\Rawfiles	APN005	Cal_01	0.000	1.000	Tray01:2	30.00	0.000	0.000
5	Standard	APN_005	\Rawfiles	APN006	Cal_01	0.000	1.000	Tray01:2	30.00	0.000	0.000
6	Standard	APN_006	\Rawfiles	APN007	Cal_02	0.000	1.000	Tray01:3	30.00	0.000	0.000
7	Standard	APN_007	\Rawfiles	APN008	Cal_02	0.000	1.000	Tray01:3	30.00	0.000	0.000
8	Standard	APN_008	\Rawfiles	APN009	Cal_02	0.000	1.000	Tray01:3	30.00	0.000	0.000
9	Standard	APN_009	\Rawfiles	APN010	Cal_02	0.000	1.000	Tray01:3	30.00	0.000	0.000
10	Standard	APN_010	\Rawfiles	APN011	Cal_02	0.000	1.000	Tray01:3	30.00	0.000	0.000
11	Standard	APN_011	\Rawfiles	APN012	Cal_03	0.000	1.000	Tray01:4	30.00	0.000	0.000
12	Standard	APN_012	\Rawfiles	APN013	Cal_03	0.000	1.000	Tray01:4	30.00	0.000	0.000
13	Standard	APN_013	\Rawfiles	APN014	Cal_03	0.000	1.000	Tray01:4	30.00	0.000	0.000

- Inspect the processing sequence and ensure that each raw data file in the three-drugs data set associates properly with a standard calibration level, QC level, unknown sample, or a blank.

Ensure that the processing sequence for the three-drugs example appears correct and proceed to [“Exercise 6: Processing the Raw Data Files and Reviewing the Analytical Results.”](#)

Saving the Workbook and Exiting the LCquan Application

If you want to take a break, you can save your workbook and open it again when you are ready to continue. For instructions about saving a workbook, see [“Saving the Workbook and Exiting the LCquan Application”](#) on page 59.

Exercise 6: Processing the Raw Data Files and Reviewing the Analytical Results

This exercise shows how to process the raw data by using the processing method and the processing sequence that you created in the previous exercises. It also explains how to review the results for each component in each raw data file and how to produce reports.

- [Processing the Raw Data Files](#)
- [Reviewing the Calibration Standards](#)
- [Reviewing the Unknowns](#)

1 Tutorial Exercises

Exercise 6: Processing the Raw Data Files and Reviewing the Analytical Results

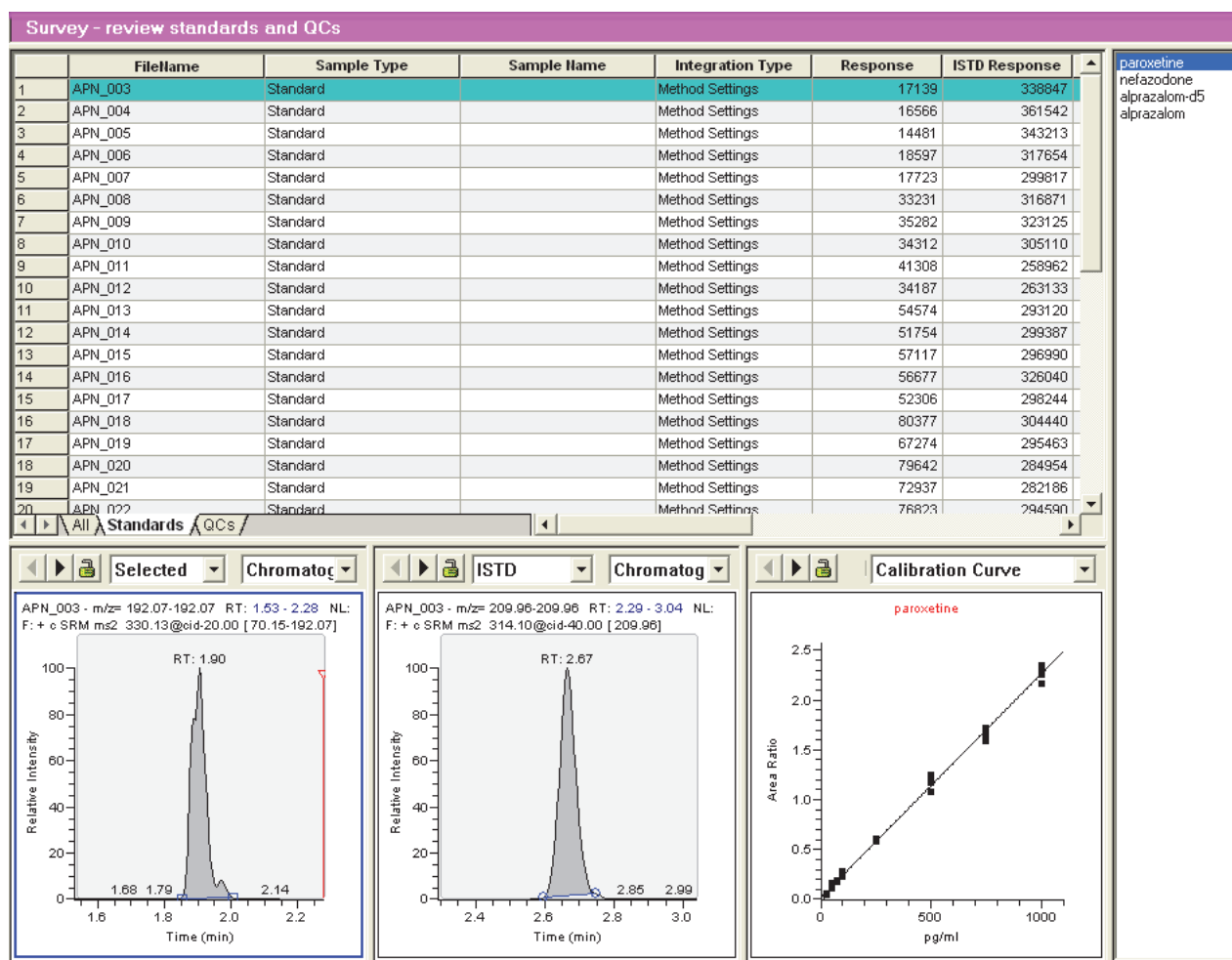
Processing the Raw Data Files

The LCQuan application processes the raw data files in the processing sequence when you open the Survey page, the Review All Results page, or the Review Reports page of the Quantitate view. The LCQuan application uses the peak detection and integration parameters in the processing method to detect and integrate peaks for all components. It then uses the calibration parameters in the processing method to build a calibration curve that it uses to quantitate the QCs and unknowns.

Although you can review the results in any order, in this example you first review the calibration standards.

Click the **Survey** button to process the raw data files and open the Survey page. In the following example (see [Figure 58](#)), the Survey page shows the results for paroxetine.

Figure 58. Paroxetine results on the Survey page



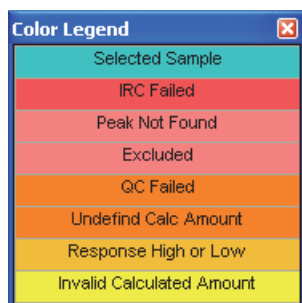
The Survey page includes (clockwise from the top) the Result List, the Component List, the Calibration Curve pane, the ISTD Chromatogram pane, and the Chromatogram pane.

The Result List contains the peak integration results. It has three display tabs: All, Standards, and QCs. Clicking the All tab displays the results for all of the standards and QCs.



The Result List also uses a system of color shading to indicate the status of individual samples (see Figure 59).

Figure 59. Color coding legend



Reviewing the Calibration Standards

❖ To review the calibration standards

1. In the Component List on the right side of the Survey page, select the **Paroxetine** target compound.

The LCquan application automatically updates the Result List, the Chromatogram and ISTD Chromatogram panes, and the Calibration Curve pane.

2. Ensure that the calibration standards are displayed in the Result List.

If necessary, click the **Standards** tab at the bottom of the Result List to display the calibration standards.

3. In the Calibration Curve pane, inspect the calibration curve at all concentrations.

When you process your own data, evaluate the calibration curve according to the criteria used in your laboratory.

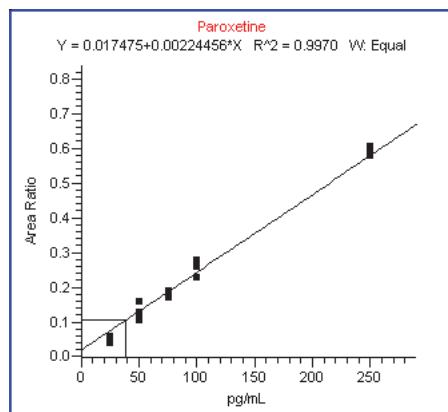
4. Zoom in to inspect the calibration curve at low concentrations, as follows:
 - a. Drag the cursor diagonally across the low concentration data points in the calibration curve.
 - b. Release the mouse button.

1 Tutorial Exercises

Exercise 6: Processing the Raw Data Files and Reviewing the Analytical Results

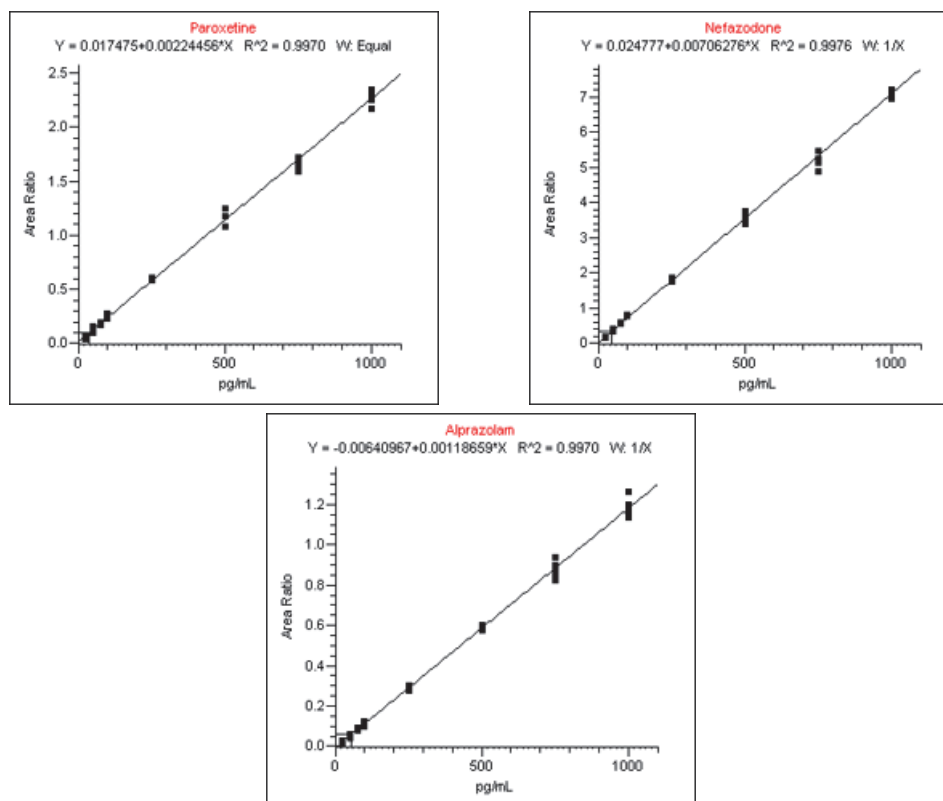
The LCQuan application zooms in on the selected data points. The Calibration Curve in [Figure 60](#) shows only the low-concentration data points.

Figure 60. Calibration curve showing only low-concentration data points



[Figure 61](#) shows the calibration curves for paroxetine, nefazodone, and alprazolam.

Figure 61. Calibration curves for the three-drugs example



5. Reset the scaling of the x and y axes of the calibration curve to full scale by clicking the **Reset Scaling**  button at the top of the page.

Tip You can change the calibration curve parameters, including curve type and level information, by using the Calibration Settings dialog box. To open the Calibration Settings dialog box, right-click the Calibration Curve pane to display a shortcut menu and choose **Calibration Settings**.

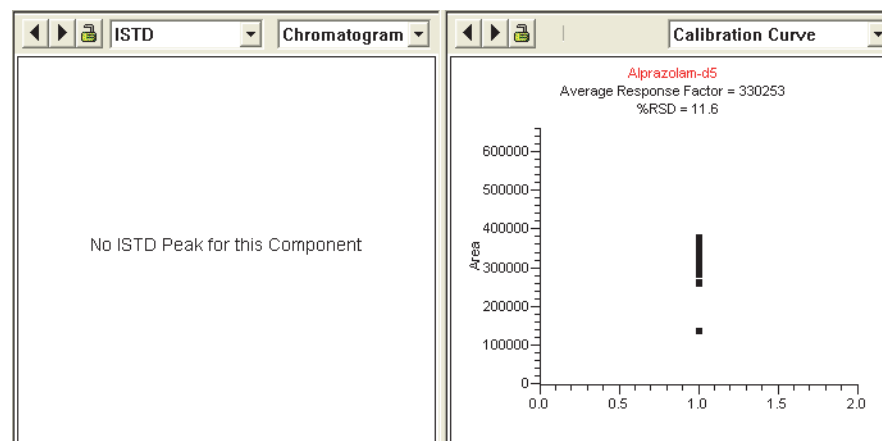
6. Inspect the Result List:
 - a. Check the entries in the Result List for peak detection and integration problems. (Use the scroll bar at the bottom of the list to see all of the entries.)
 - b. Ensure that the data files correspond to the correct levels and sample types.
7. Select the first data file by clicking the first row of the Result List.
8. Inspect the component and internal standard peaks, as follows:
 - a. Inspect the component peak in the Chromatogram pane:
 - i. Ensure that the LCQuan application found the peak.
LCQuan shades the peak gray and marks the starting and ending points of the peak with integration markers.
 - ii. Ensure that the shaded area accurately represents the contribution of the component to the chromatogram.
 - b. Inspect the internal standard peak in the ISTD Chromatogram pane:
 - i. Ensure that the LCQuan application found the peak.
LCQuan shades the peak gray and marks the starting and ending points of the peak with integration markers.
 - ii. Ensure that the shaded area accurately represents the contribution of the internal standard to the chromatogram.
9. Inspect the remaining files in the Result List:
 - a. Select the next data file by clicking on the next row in the Result List.
 - b. Repeat step 8 for each file.
10. Inspect the ion ratio chromatograms for nefazodone and paroxetine:
 - a. In the Components list, select the component.
 - b. In any of the bottom panes, select **IRC Chromatogram** from the list.
11. Review the internal standard results as you did for the target compound.
 - a. In the Component List, select the internal standard, **Alprazolam-d5**.

The LCQuan application automatically updates the Result List and the Chromatogram pane. LCQuan does not display a chromatogram in the ISTD Chromatogram pane for this component.

1 Tutorial Exercises

Exercise 6: Processing the Raw Data Files and Reviewing the Analytical Results

The LCQuan application displays the response for each internal standard sample in the Calibration Curve pane.



- b. Ensure that the data files correspond to the correct levels and sample types.
- c. Inspect the chromatogram in the Chromatogram pane for each file:
 - Ensure that the LCQuan application found the peak.
 - Ensure that the shaded area accurately represents the contribution of the component to the chromatogram.

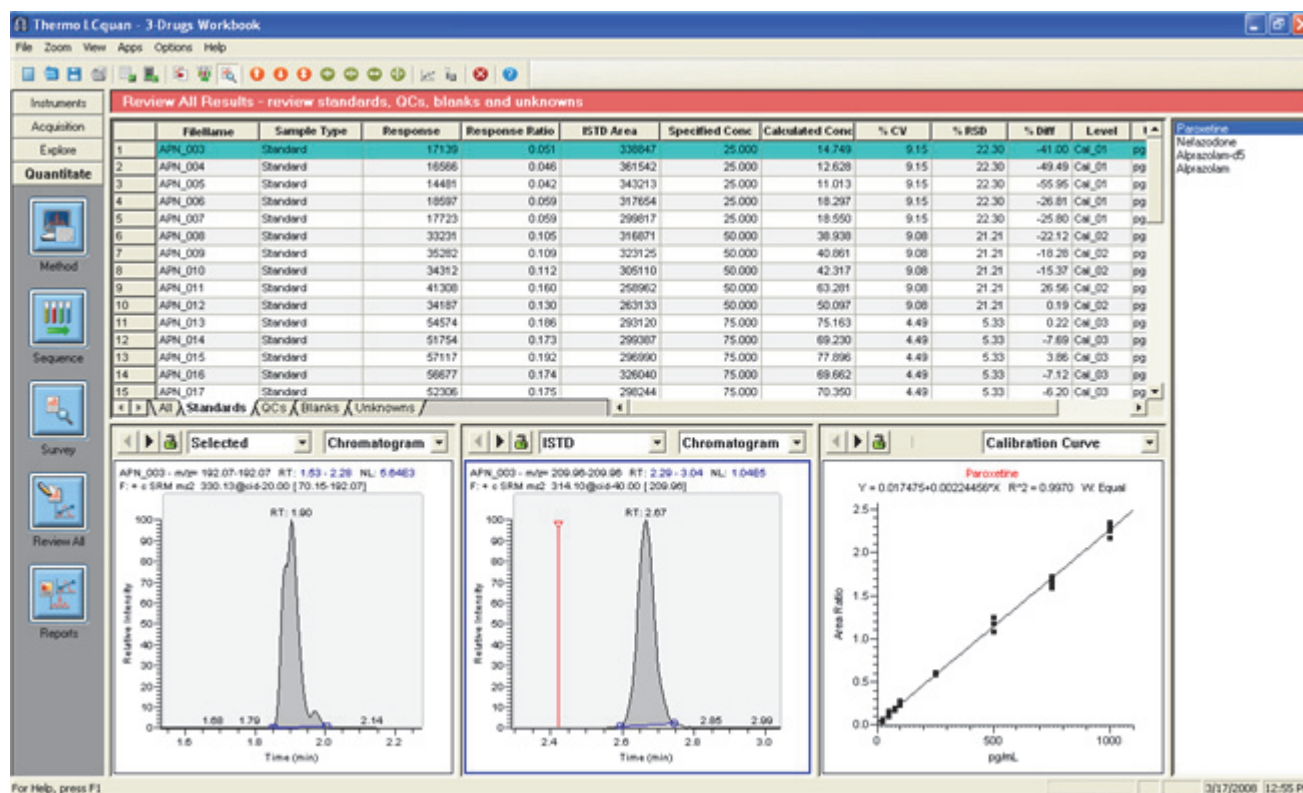
Reviewing the Unknowns

Use the Review All Results page on the Quantitate view to review and rework samples of all types, including unknowns and blanks. For calibration standards and QCs, the Review All Results page functions exactly as the Survey page does.

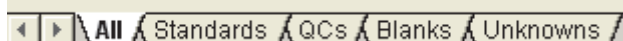
To process the raw data files and open the Review All Results page, click the **Review All** button in the navigation pane.

Figure 62 shows the results for the three-drugs data.

Figure 62. Review All Results page in the Quantitate view



The Review All Results page has five display tabs: All, Standards, QCs, Blanks, and Unknowns. Clicking the All tab displays the results for every sample in the sequence.



❖ To review and rework the results for the unknowns

1. Click the **Unknowns** tab at the bottom of the Result List to display the peak integration results for the unknown samples.
2. In the Component List, select the internal standard, **Alprazolam-d5**.
3. Inspect each file in the Result List and check for peak detection and integration problems.
4. Inspect the chromatogram in the Chromatogram pane for each file, as follows:
 - a. Ensure that the LCquan application found the peak.
 - b. Ensure that the shaded area accurately represents the contribution of the component to the chromatogram.

If you want to modify the peak detection and integration settings, return to “[Entering Component Identification Parameters and Integrating the Peaks](#)” on [page 32](#).

1 Tutorial Exercises

Exercise 6: Processing the Raw Data Files and Reviewing the Analytical Results

5. In the Component List, select a target compound.
6. Repeat steps 3 and 4 for each target compound.

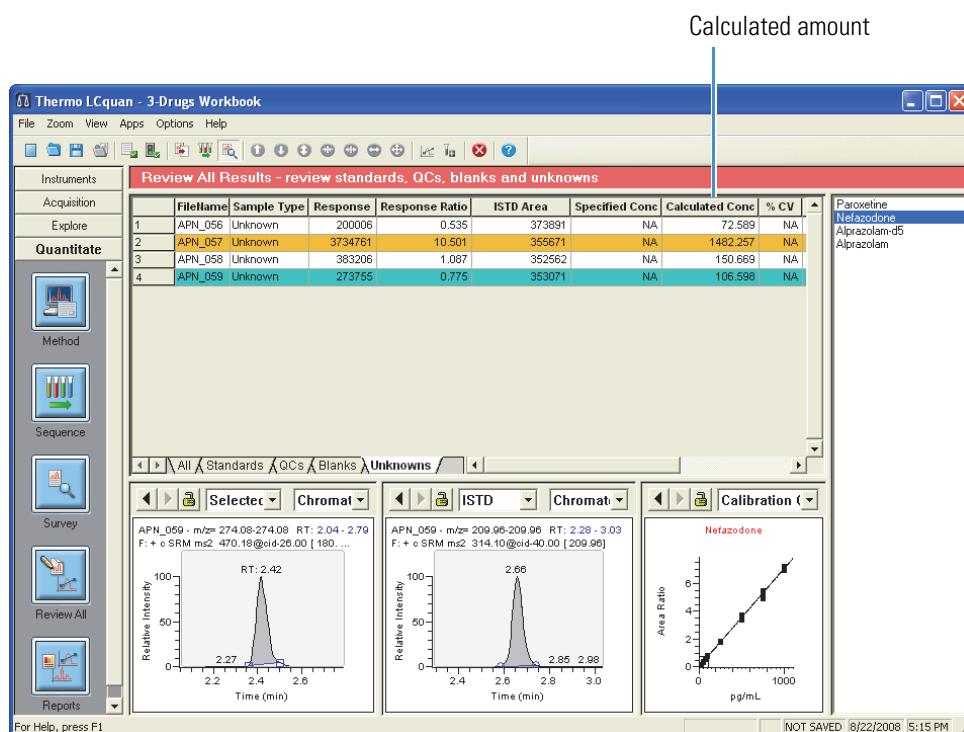
Note For the file APN_057, target compound **nefazodone**, the row in the Result List is shaded in pale orange and the Peak Status column displays **Response High**. This means that the area ratio for the sample is greater than the area ratio for any of the calibration standards. As a result, the LCquan application had to extrapolate the calibration curve to determine the amount.

7. After confirming that the LCquan application has detected and integrated the peaks properly, inspect the Result List again.

The Calculated Amount column of the list displays the calculated amount of drugs in each unknown sample.

Figure 63 shows the results for the nefazodone unknowns.

Figure 63. Results for nefazodone unknowns on the Review All Results page



8. Go to “Exercise 7: Generating a Custom Report.”

Saving the Workbook and Exiting the LCquan Application

If you want to take a break, you can save your workbook and open it again when you are ready to continue. For instructions about saving a workbook, see “Saving the Workbook and Exiting the LCquan Application” on page 59.

Exercise 7: Generating a Custom Report

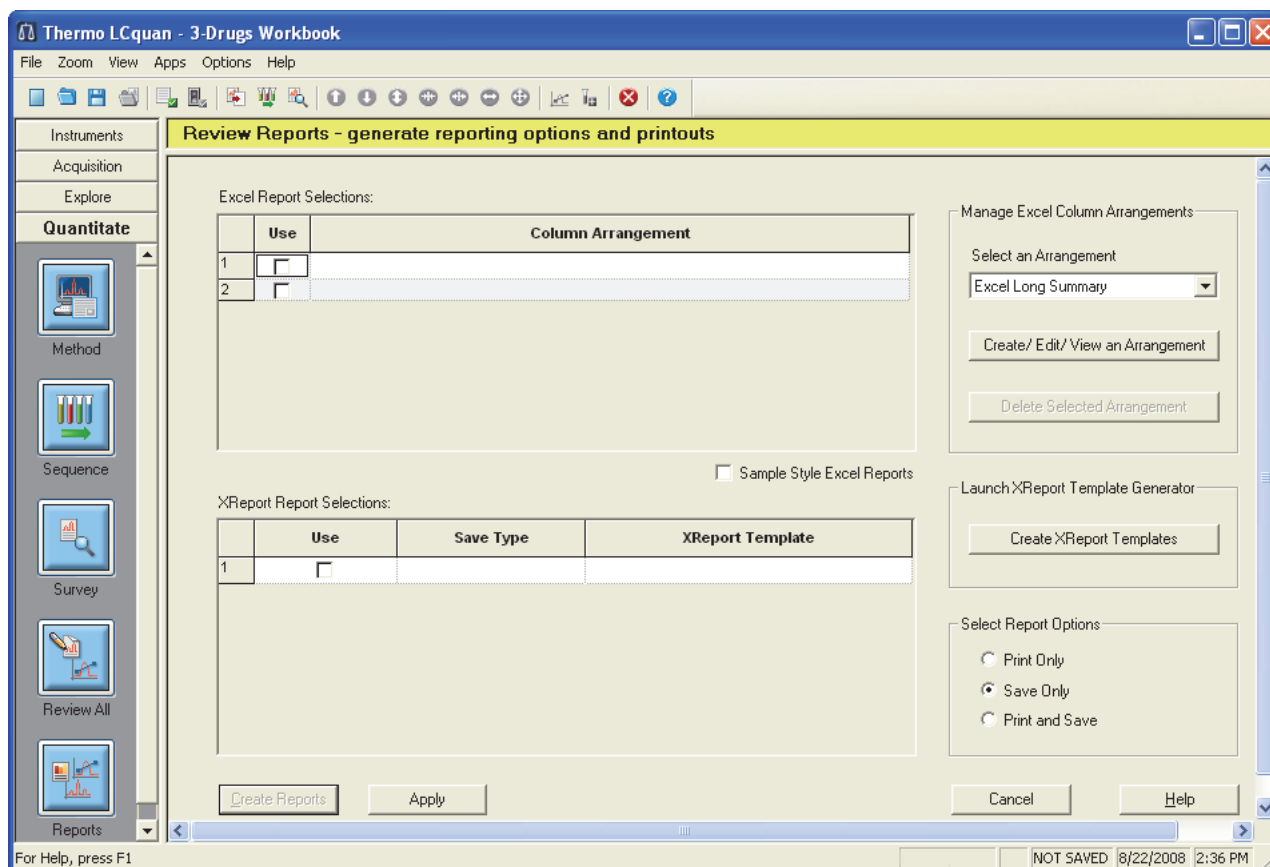
This exercise describes how to create a custom Microsoft Excel™ report that describes the quantitative analysis results for your data.

❖ To generate a custom report

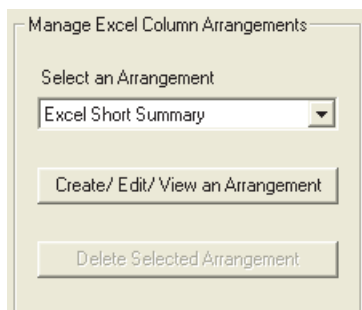
1. Click the **Reports** icon in the navigation pane.

The Review Reports page opens (see [Figure 64](#)).

Figure 64. Review Reports page in the Quantitate view

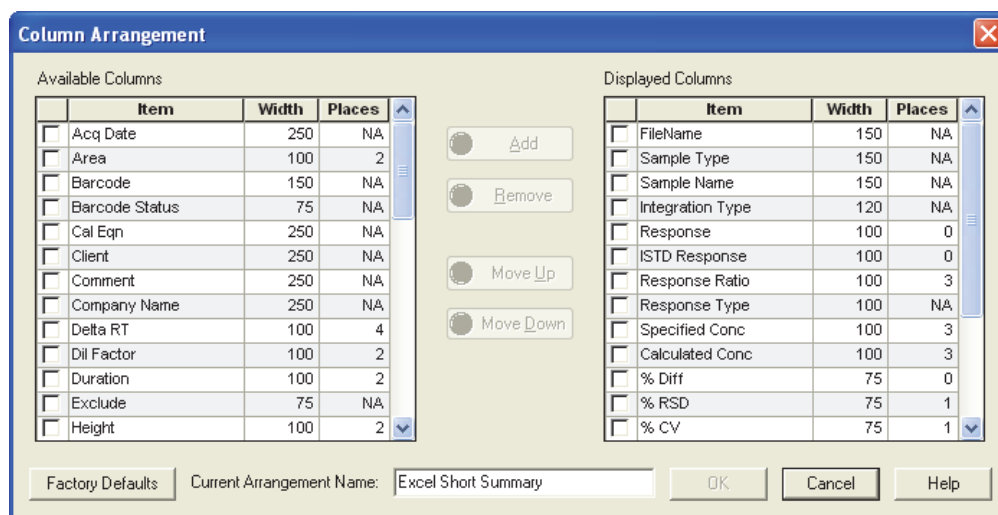


2. In the Manage Excel Column Arrangements area (see [Figure 65](#)), select **Excel Short Summary** from the list.

Figure 65. Mange Excel Column Arrangements area

3. Click **Create/Edit/View an Arrangement**.

The Column Arrangement dialog box opens (see [Figure 66](#)).

Figure 66. Column Arrangement dialog box

The Displayed Columns list shows all the columns that are included in the Excel Short Summary report. The Available Columns list shows (in alphabetical order) all the available columns that are not included in this report style.

You cannot modify the standard Excel Short Summary report, but you are going to use it as a template to create your custom report.

4. Change the Current Arrangement Name to **CustomSummary**.

Notice that the OK button is enabled so that you can now save your changes to this report style.

5. In the Available Columns list, click the box before Acq Date and click **Add**.

The Acquisition Date column is added to the columns that will be displayed in your report.

6. In the Displayed Columns list, click the boxes before any of the columns you do not want in your report and click **Remove**.

7. Use the **Move Up** and **Move Down** buttons to reorder the displayed columns any way you like.
8. When you have finished editing the columns for the new arrangement, click **OK**.
9. In the Excel Report Selections table (see [Figure 67](#)), select the **Use** check box in an empty row.

Figure 67. Excel Report Selections area

	Use	Column Arrangement
1	☒	CustomSummary
2	☐	Quan Result Grid Excel Long Summary Excel Short Summary CustomSummary

☐ Sample Style Excel Reports

10. Click the **Column Arrangement** column and select **CustomSummary** from the list.
11. To create sample style Excel reports, select **Sample Style Excel Reports**.
Sample Style Excel Reports have information about one sample per page.
12. In the Select Report Options area (see [Figure 68](#)), select the **Save Only** option.

Figure 68. Select Report Options area

Select Report Options

☐ Print Only
☒ Save Only
☐ Print and Save

13. Click **Create Reports**.
14. When asked if you want to save your workbook, click **Yes**.

The application writes the generated Excel (.xls) report to the Exports folder:

...\\Xcalibur\\QuanRoot\\Example Study\\3-Drugs Workbook\\Exports

Saving the Workbook and Exiting the LCquan Application

You can save the workbook at any step in the LCquan procedure. You can then exit the LCquan application and reenter at the same place at a later time. When you reopen a workbook, LCquan displays the page and view that were open when you exited the workbook.

1 Tutorial Exercises

Saving the Workbook and Exiting the LCQuan Application

❖ To save the workbook and exit the LCQuan application

1. Choose **File > Save** to save the current workbook.

If you want to save the workbook with a different study or workbook name, choose **File > Save As**. The Save As Wizard guides you through the procedure.

2. Choose **File > Exit** to exit the LCQuan application.

Note During data acquisition, you cannot exit the application without saving the workbook.

Index

A

acquisition sequence, editing [14](#)

C

calibration

curve parameters, specifying [30](#)

levels, assigning [18](#)

standards, reviewing [51](#)

chromatogram peaks, relating to components [22](#)

components, naming [18](#)

contacting us [vi](#)

creating a new study and workbook [2](#)

creating a processing method [27](#)

custom report, generating [57](#)

D

designing a custom report [57](#)

downloading documents [vi](#)

E

editing, acquisition sequence [14](#)

exercises

create acquisition sequence [7](#)

create new study and workbook [2](#)

create processing method [27](#)

create processing sequence [42](#)

explore data [19](#)

generate custom report [57](#)

process raw files and review results [49](#)

Explore

results, description [23](#)

Review page [23](#)

G

generating a custom report [57](#)

I

identification settings, specifying [32](#)

Import Data page [7](#)

Import from Existing Workbook parameter [6](#)

Import Legacy Files parameter [6](#)

internal standard, identifying [29](#)

N

new study (and workbook), creating [2](#)

P

peak detection, description [21](#)

processing method, creating [27](#)

processing sequence, reviewing [48](#)

R

raw data files

associating with sample types [43](#)

opening [20](#)

processing [50](#)

Rawfiles folder, browser [5](#)

S

scan filters [21](#)

T

Thermo Fisher website, user documents [vi](#)

U

unknowns, reviewing [54](#)

W

welcome page [3](#)

Wizards

New Acquisition Sequence [8](#)

Index: W

New Method [27](#)
New Processing Sequence [42](#)
New Study or Workbook [4](#)
workbook, saving [60](#)