



Agilent 6530 LC/Q-TOF

Introduction Workbook



Notices

Document Identification

G1960-90153

January 2026

Copyright

© Agilent Technologies, Inc.

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Agilent Technologies, Inc.
5301 Stevens Creek Blvd.
Santa Clara,
CA
95051
www.agilent.com

Software Revision

This workbook is valid for MassHunter 12.2, until superseded.

Warranty

The material contained in this document is provided “as is,” and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

Safety Notices

CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

Contents

1	Introduction	5
	About this Training Workbook	6
	How to use this Workbook	7
	Before you begin	7
	Additional Resources	8
	Where to find more information	8
2	Hardware	9
	Overview	10
	Front view	10
	Side view	11
	Back view	12
	Basic Components	13
	Ionization Source	13
	Calibrant Delivery System	15
	Nebulizer	16
	Rough Pump	17
	Waste Bottle from LC Pump or Mass Spectrometer	18
3	Software Basics	19
	Software Start-Up	20
	User Interface and General Navigation	21
	Close Connection	22
	Creating and Configuring Projects	23
	Instruments	24
	Create an Instrument	25
	Launch an Instrument	26
	Create an instrument shortcut	27
4	Tuning	28
	What is tuning in LC/MS?	29
	What is the difference between Tunes?	29
	Mass Calibration	31

	Calibrate the LC/Q-TOF (checktune)	32
	Review a Tune Report	33
	Schedule a Tune	34
	Stop checktune	35
5	Using Methods	36
	The Method Editor Window	37
	Set the Active Instrument Mode	38
	Set Up and Run an Acquisition Method	39
	Load the default method	39
	Set Up a Scan Method	42
	Run Methods	46
	Review the Data using MassHunter Qualitative Analysis	49
6	Run a Worklist	52
	Create and edit a worklist	53
	Add Multiple Samples	54
	Change Result Window Layouts	56
	Use layout icons on the toolbar	56
	Change the panes in the Compound Information window for Cal-L4	57
	Save the default layout without the calibration curve	58
	Load the newly created layout	59
	Recreate (do not restore) the default layout	61
7	Maintenance	62
	Register on Agilent SubscribeNet (New Account Registration)	63
	Perform Back Up and Platform Best Practices	64



1 Introduction

About this Training Workbook	6
How to use this Workbook	7
Before you begin	7
Additional Resources	8
Where to find more information	8

About this Training Workbook

This workbook provides instructions for the Agilent 6530 LC/Q-TOF system running MassHunter Acquisition 12.2 or higher.

For more information on the software and detailed instructions on the workflow not covered in this workbook, see the Online Help.

This workbook is your introductory guide for the set-up and execution of basic procedures with the LC/Q-TOF and method development workflow. This workbook is divided into chapters, each building upon the last, so we recommend that each chapter is completed in succession. During each chapter, lessons are guided by an Agilent-certified service professional.

By completing this learning event, you will have an introductory level of experience in the use of an Agilent 6530 Q-TOF LC/MS system.

This introduction covers:

- Reviewing hardware components and software procedures.
- Performing a checktune.
- Acquiring and analyzing a sample.
- Performing routine maintenance.

How to use this Workbook

This learning experience introduces basic concepts in a learning-by-doing, guided manner. Each chapter uses step-by-step instructions.

Exercises to be completed are marked like this:



Exercise Name

Exercise Instructions

Task steps look like this:

1. Tasks or items needed to complete tasks look like this.

If you are expected to enter any information or if something is important, it is set in italicized type like this:

Type *Blank One* in the field.

If you are expected to press a key on the keyboard or button on the software screen, the key is displayed in bold like this:

Press **Enter**.

Before you begin

This introduction workbook is recommended for all participating end users.



- Download the *Agilent 6530 Q-TOF LC/MS System User Guide* by scanning the code or navigating to <https://aglt.co/LCMSUserDocs>.

Introduction

Where to find more information

- Use the *Agilent 6530 Q-TOF LC/MS System Introduction Workbook* and Introduction Checklist with your Agilent-certified service professional and keep for future reference.

Additional Resources

User Documentation



Data analysis and library management documentation can be found by scanning the code or navigating to <https://aglt.com/DALibMgmtDocs>.



Instrument documentation, step-by-step videos, and more can be found by scanning the code or navigating to <https://aglt.co/LCMSUserDocs>.

Agilent Q-TOF LC/MS Supplies



Use this quick reference list to keep your shelves stocked by navigating to <https://aglt.co/LCQTOFSupplies>.

Where to find more information



Agilent Community

To get answers to your questions, join over 10,000 users in the Agilent Community. Review curated support materials organized by platform technology. Ask questions to industry colleagues and collaborators. Get notifications on latest videos, documents, tools, and webinars relevant to your work.

<https://community.agilent.com/>



2 Hardware

Overview	10
Front view	10
Side view	11
Back view	12
Basic Components	13
Ionization Source	13
Calibrant Delivery System	15
Nebulizer	16
Rough Pump	17
Waste Bottle from LC Pump or Mass Spectrometer	18

Overview

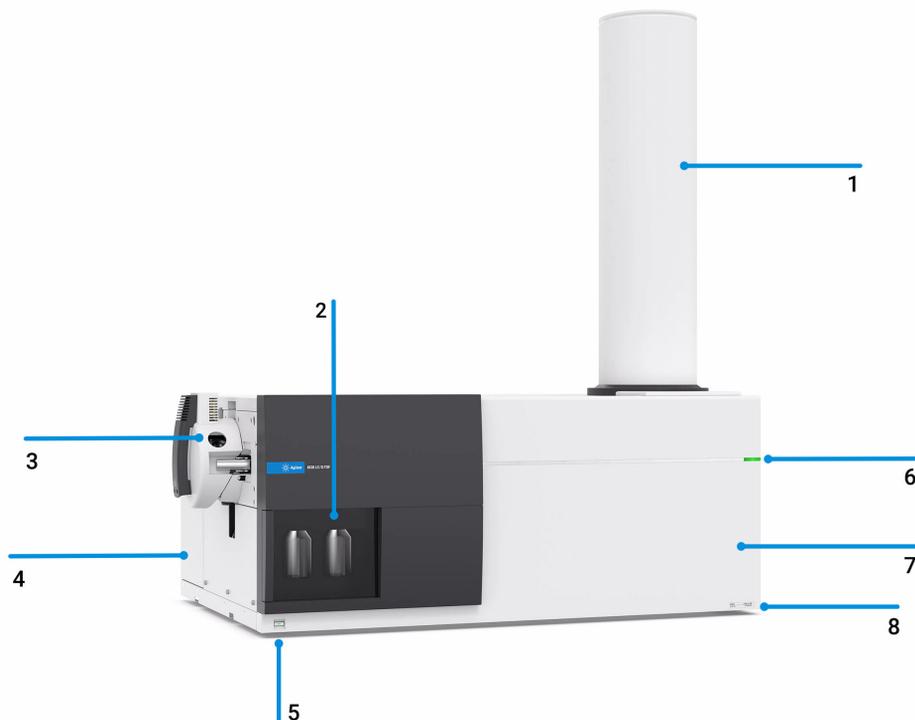
In this section, you will identify basic hardware components and their locations for the Agilent 6530 LC/Q-TOF system.



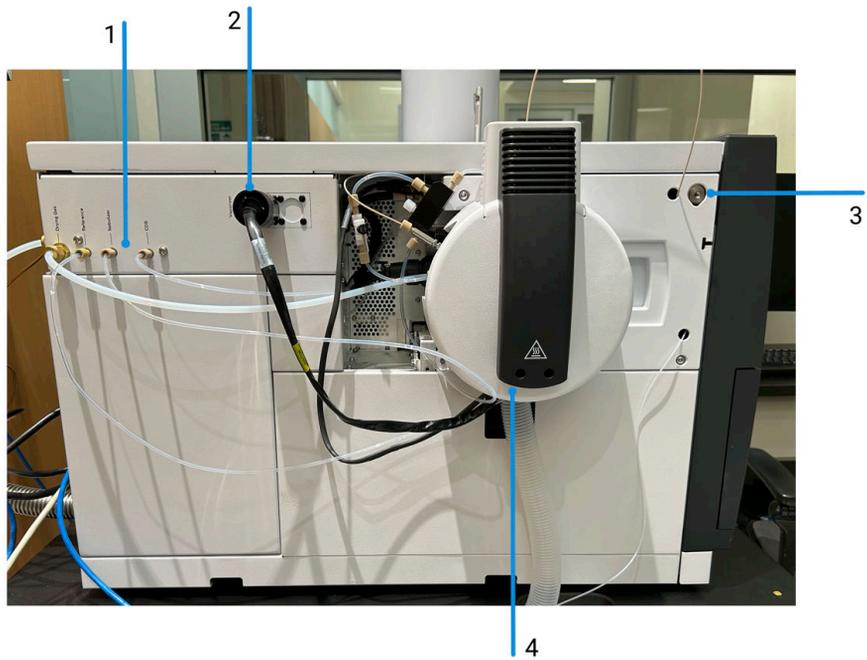
Fill in the Blank

Fill in the Blank: Work with your Agilent Service Engineer and/or use the *Agilent 6530 LC/Q-TOF System User Guide* to label the flagged components below for your installed instrument(s).

Front view



Side view



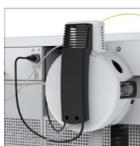
Back view



Basic Components

Ionization Source

Agilent liquid chromatography/mass spectrometry (LC/MS) ion sources enable analysis of a wide range of samples quickly and accurately. These easily interchangeable ion sources for Agilent LC/MS systems include the:



- Agilent Jet Stream (Dual AJS) source



- Electrospray Ionization (Dual ESI) source



- Atmospheric Pressure Chemical Ionization (APCI) source



- Multimode Ionization source (MMI)

Hardware

Basic Components



- Agilent nanospray ESI source



Hardware Introduction

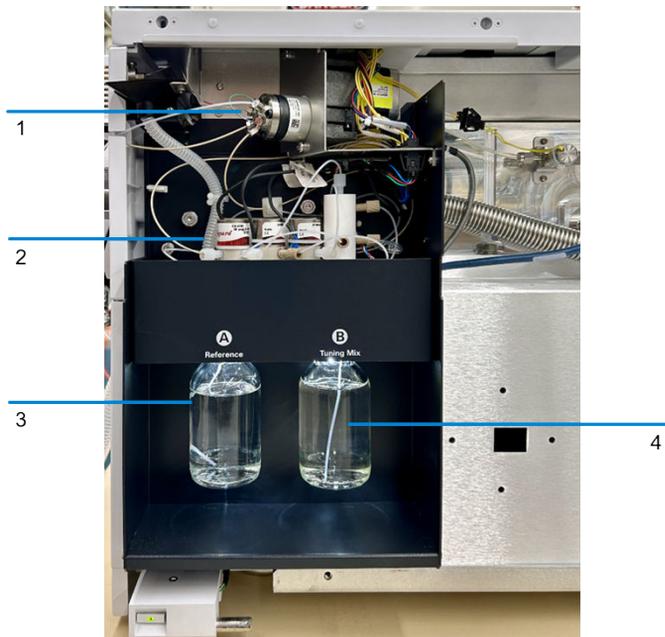
1. List the ionization source in use:

2. It is reviewed on page ____ of the user guide and includes the following parts:

3. List the name and part number of the proper tune solution for this system:

Calibrant Delivery System

The calibrant delivery system (CDS) introduces calibration solution for automated mass calibration of the mass spectrometer, to ensure that the mass accuracy of the system is maintained throughout batch acquisition.



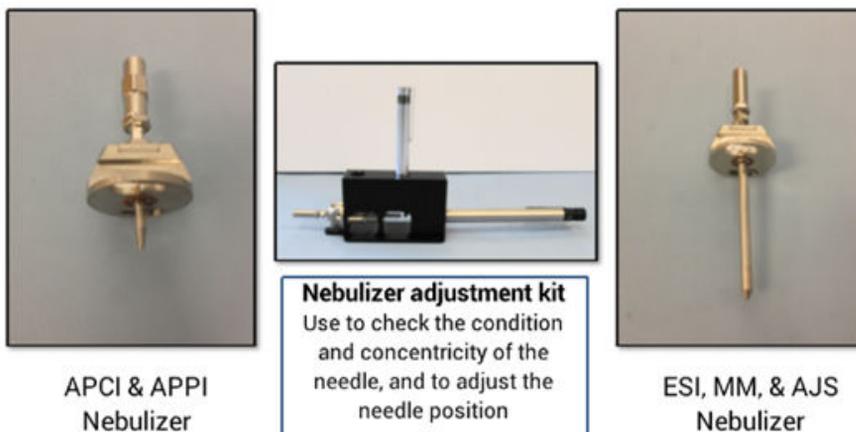
Hardware Introduction

1. Practice removing and attaching the calibrant bottle and the internal reference mass bottle.
2. Are the calibrant bottle and the internal reference mass bottle interchangeable? Why?

3. How often is the calibrant bottle checked and refilled?

Nebulizer

A nebulizer is a device for producing a fine mist of charged droplets that converts a liquid sample into an aerosol for introduction into the vacuum system.



View the Needle

- Find your nebulizer type per the user guide or the document that comes with the kit. List the part number below:

Rough Pump

MS40+



Locate the Oil Sight

Using the user guide, fill out the following information:

1. The oil level should be _____ the marks for Max and Min.
2. Check that the pump oil is _____ and the color is _____ than amber.
3. If the pump oil is _____ or full of _____, replace it.

Waste Bottle from LC Pump or Mass Spectrometer



Waste Bottle

1. At what level should you empty the waste bottle?

2. Does the waste bottle have a waste line connected? Why is this important to keep in place?



3 Software Basics

Software Start-Up	20
User Interface and General Navigation	21
Close Connection	22
Creating and Configuring Projects	23
Instruments	24
Create an Instrument	25
Launch an Instrument	26
Create an instrument shortcut	27

Software Basics

Software Start-Up

The MassHunter Control Panel is the administrative and management center for MassHunter Data Acquisition software. Control Panel provides:

- Full instrument status information of your entire laboratory
- Central configuration and administration of users, instruments, and security settings
- Full system documentation and built-in reports

You will review:

- Starting the software
- Navigation overview
- Closing the connections
- Creating projects
- Creating and configuring instruments
- Launching instruments
- Offline method editor
- Creating shortcuts

Software Start-Up



- From the desktop, double-click the **Control Panel** icon.



The navigation pane opens by default and can be minimized or expanded based on your preference.

User Interface and General Navigation

Ribbon

Panes
Instruments - Controls specific instruments.
Projects - Create paths to save project data.
Administration - Add and remove configuration

Main Window

Adjusting the Instruments pane:

- To minimize the pane, click <<. When minimized, the tab currently selected is displayed vertically.
- To expand the pane, click >>.
- Items can be dragged and dropped in the Instruments and Projects pane. The existing privileges of the instrument or project are not retained when moving. The user must have the proper privileges to perform this function.

Close Connection

Use the Close Connection function to sever the connection between the instrument and the configured Instrument Controller (AIC or Workstation).

1. Click **Instruments**.
2. Select the instrument to close.
3. Click **Close Connection**.

The screenshot shows the 'LC/Q-TOF - Control Panel' window. The top management bar contains several icons, with 'Close Connection' highlighted by a yellow box. The left sidebar has 'Instruments' selected, with an orange arrow pointing to it. Below it, the 'LC/Q-TOF' instrument is selected, with another orange arrow pointing to it. The main panel displays the instrument's status as 'Idle' and an activity log table.

Date/Time	User	Description
2023-11-29 08:54:22-08:00	SYSTEM (SYSTEM)	Instrument "Instruments\Revident LC/Q-TOF" was changed
2023-11-29 08:09:38-08:00	SYSTEM (SYSTEM)	Acquisition Engine: Method 'C:\Projects\Learning Products\Method
2023-11-29 08:09:18-08:00	SYSTEM (SYSTEM)	Acq Console: User 'SYSTEM' has logged in.
2023-11-29 08:09:15-08:00	SYSTEM (SYSTEM)	Engine Launcher: System Engines started.
2023-11-29 08:08:13-08:00	SYSTEM (SYSTEM)	Configuration of instrument "Instruments\DESKTOP-655PFQM" was
2023-11-29 08:08:13-08:00	SYSTEM (SYSTEM)	Instrument Configuration: User 'SYSTEM (SYSTEM)' has logged out.
2023-11-29 08:08:13-08:00	SYSTEM (SYSTEM)	Instrument "Instruments\DESKTOP-655PFQM" was configured
2023-11-29 08:07:48-08:00	SYSTEM (SYSTEM)	Instrument Configuration: User 'SYSTEM (SYSTEM)' has logged in.
2023-11-29 08:06:17-08:00	SYSTEM (SYSTEM)	Instrument Configuration: User 'SYSTEM (SYSTEM)' has logged out.

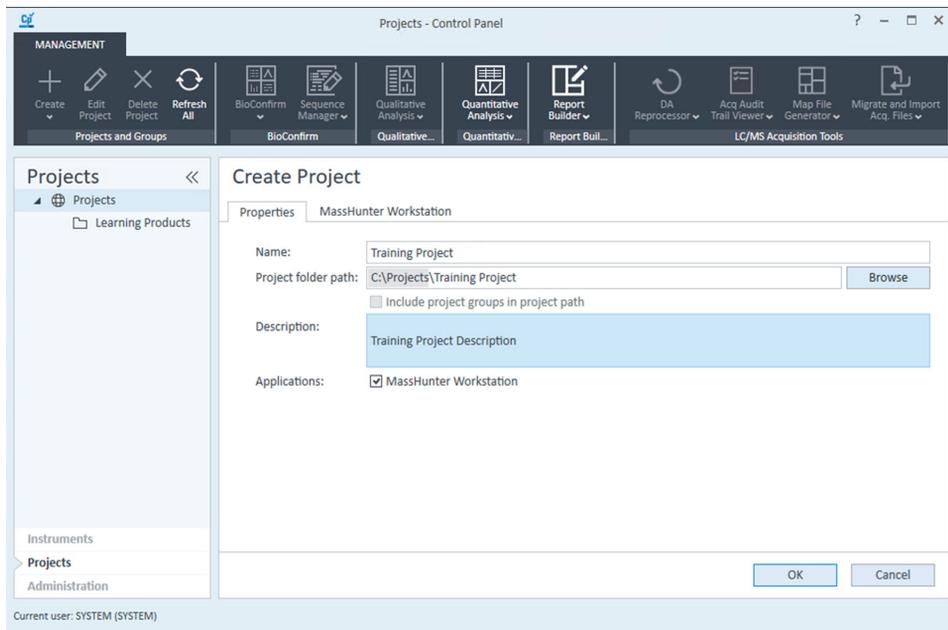
Current user: SYSTEM (SYSTEM)

Creating and Configuring Projects



Creating and Configuring Projects

1. Click **Projects** and select **Projects**.
2. In the **Name** text box, type *Training Project*.
3. In the Project folder path text box, leave the default folder path.
4. In the Description text box, type a description of the project, for this example *Training Project Description*.
5. Click the **MassHunter Workstation** tab and review the available options. Do not change the defaults.
6. Click **OK**.



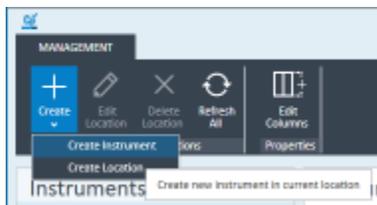
Instruments

Use the Control Panel to connect and control the instruments you want to use with the software.

Create an Instrument



1. Click **Instruments** and select any location.
2. Click **Create > Create Instrument**.

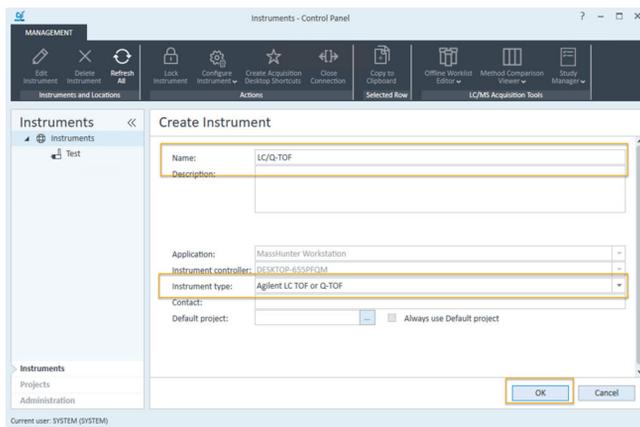


3. Enter the required data in the Create Instrument pane.
 - a. Name: 6530 LC/Q-TOF
 - b. Instrument Type: **Agilent LC TOF or Q-TOF**

NOTE

Do not select a default project. You will be prompted to select a project when you launch the instrument.

4. Click **OK**.
5. Click ... to select the **TrainingProject** project from the Select Project dialog.
6. Click **OK**. The instrument is displayed in the navigation pane.



Launch an Instrument

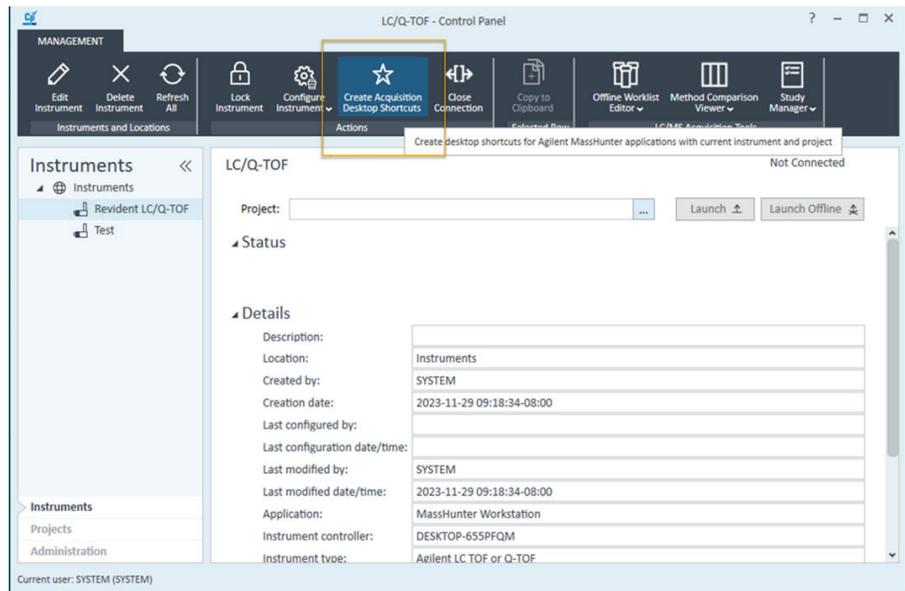
Once you have added an instrument, launch the instrument to begin acquisition from the instrument table or the instrument details page, or launch an instrument directly from your desktop shortcut.

1. Click **Instruments** and select an instrument from the left panel.
2. In the Instruments window, click **Launch**.

The screenshot shows the 'LC/Q-TOF - Control Panel' window. The top toolbar includes icons for 'Edit Instrument', 'Delete Instrument', 'Refresh All', 'Lock Instrument', 'Configure Instrument', 'Create Acquisition Desktop Shortcuts', 'Close Connection', 'Copy to Clipboard', 'Offline Worklist Editor', 'Method Comparison Viewer', and 'Study Manager'. The left sidebar shows a tree view with 'Instruments' selected, and 'LC/Q-TOF' is highlighted. The main content area displays the instrument details for 'LC/Q-TOF', including a 'Project' field, a 'Launch' button (highlighted with a yellow box), and a 'Launch Offline' button. Below this is an 'Activity Log (last 7 days)' table with one entry: '2023-11-29 09:18:34-08:00' by 'SYSTEM (SYSTEM)' with the description 'Instrument "Instruments\Revident LC/Q-TOF" was added'. The status at the top right is 'Not Connected'. The current user is 'SYSTEM (SYSTEM)'.

Create an instrument shortcut

1. In the Control Panel, click **Instruments** and select an instrument from the left panel. Verify that the correct Project is selected.
2. Click **Create Acquisition Desktop Shortcuts** in the **Actions** group on the ribbon. Two icons are added to the desktop with the name of the instrument and whether it is online or offline.





4 Tuning

What is tuning in LC/MS?	29
What is the difference between Tunes?	29
Mass Calibration	31
Calibrate the LC/Q-TOF (checktune)	32
Review a Tune Report	33
Schedule a Tune	34
Stop checktune	35

Tuning

What is tuning in LC/MS?

When the LC/MS Q-TOF is used as a detector for the LC, a mass spectrum is associated with each data point in the LC chromatogram. To obtain high quality, accurate mass spectra, the LC/MS Q-TOF must be optimized to:

- Maximize sensitivity
- Maintain acceptable resolution
- Ensure accurate mass assignment

What is tuning in LC/MS?

Tuning is the process of adjusting LC/MS Q-TOF parameters to achieve the optimized goals listed above.

Tuning acts as a diagnostic tool to indicate the service or cleaning requirements of the spectrometer; it provides a chronicle of system performance and the matching of fragments from a known calibration compound to adjust the mass axis so it agrees with the expected mass assignments.

What is the difference between Tunes?

Mass Calibration performs a TOF mass calibration on the Active Instrument Mode, that is displayed in the Systems Settings panel. Run Mass calibration daily, at a minimum weekly. During normal operation, it take approximately 15 minutes to complete.

A checktune is run each day an analysis is performed. A checktune can be used to determine if the tuning mix ion masses are properly assigned and if the response or sensitivity of these ions is within expectations. In other words, a checktune performs a single profile scan of the tune masses and compares the peak widths and mass axes with target values to make sure they are correct before you start your acquisition. Checktune can be performed in either positive or negative ionization mode, or both. During normal operation, it take approximately 25 minutes to complete.

Tuning

What is tuning in LC/MS?

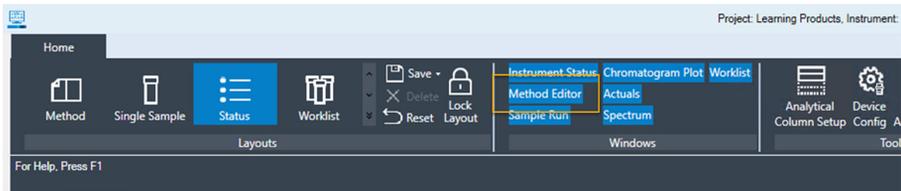
Perform an autotune monthly, after preventative maintenance, or if you find a problem with checktune. Autotune can be performed in negative and/or positive ionization.

After pumpdown, wait ~12 hours for the analyzer to reach thermal equilibrium before before tuning or operating your LC/Q-TOF system. Tune files that are created, or data that is acquired, before the LC/Q-TOF system is at thermal equilibrium may have incorrect mass assignments and other inaccuracies.

Mass Calibration

Mass calibration is run with the following ion sources: ESI, AJS ESI, and APCI.

1. In MassHunter Acquisition, click **Method Editor**.



2. Click the **MS Q-TOF** tab.



3. Click **Tune > Mass Calibration** in the left pane.
4. Click **Tune control**.
5. Select **Mass Calibration**.



6. Click **Mass calibration/Checktune**.
7. When the tune completes, review the report.
8. Click **Release tune control** in the toolbar to release control of the instrument.

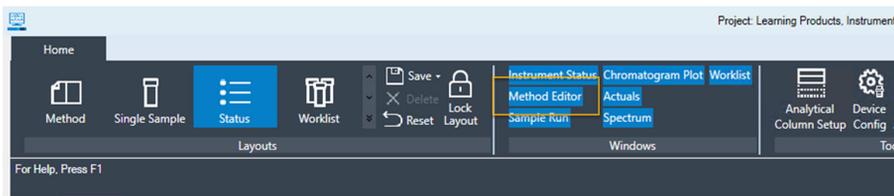
Tuning

Calibrate the LC/Q-TOF (checktune)

Calibrate the LC/Q-TOF (checktune)

A checktune is run with the following ion sources: ESI, AJS ESI, and APCI.

1. In MassHunter Acquisition, click **Method Editor**.



2. In the Method Editor window, click the **MS Q-TOF** tab.



3. Click **Tune > Mass Calibration** in the left pane.
4. Click **Tune Control**.
5. Select **Checktune**.



6. Click **Mass calibration/Checktune**.
7. When the tune completes, review the report.
8. Click **Release tune control** in the toolbar to release control of the LC/Q-TOF instrument.



Review a Tune Report

Generate a detailed tune report after you have run autotune or checktune.

1. In the Method Editor window, select the **MS Q-TOF** tab.
2. Click the **Tune > Autotune** section in the left pane.
3. Click  **Tune control**.
4. Click  **Tune Reports** in the toolbar in the Autotune section.
5. Select the tune report to view and click  **View tune reports**. If no reports are available, then run either a mass calibration, checktune, or autotune first.
6. Click  to release control of the instrument.

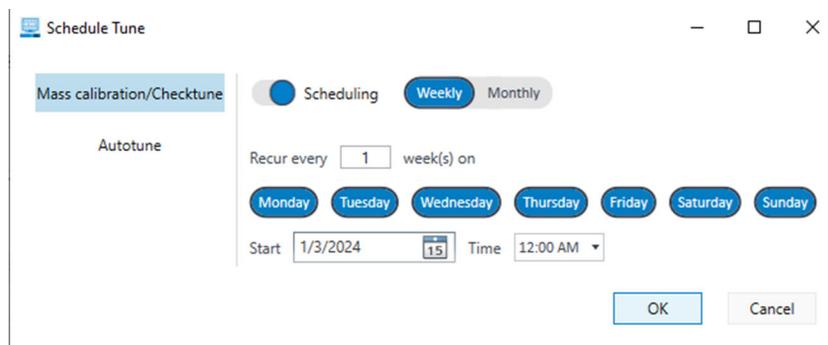
NOTE

Only the polarities that were last calibrated or tuned appear in the tune reports saved with the data files. Prior detailed tune reports can be accessed through the Autotune or Checktune section. For all other instruments, save tune reports manually before tuning in a single polarity.

Schedule a Tune

Schedule a mass calibration, checktune, or autotune so that a tune is run automatically at specified times.

1. To display the Method Editor window, click **Method Editor** in the Windows section on the Ribbon. Or, click the **Method** layout on the Ribbon.
2. In the Method Editor window, select the **MS Q-TOF** tab.
3. Select the **Tune > Autotune** section.
4. Click  **Tune control** in the toolbar to lock control of the instrument. You can't start a single sample run or a worklist when Tune has control of the instrument.
5. Click  **Schedule tune** in the toolbar.
6. Select **Mass calibration/Checktune** in the left pane. The right pane shows the information for scheduling a checktune.
7. Click the **Scheduling** slider to switch on Scheduling.
8. Select **Weekly** for this exercise.
9. Select a day of the week and a Start date and time to indicate how often to schedule the checktune.



10. Click **OK**.

Stop checktune

1. Click the **Tune > Mass Calibration** section in the left pane.
2. Click  **Tune control** in the toolbar.
3. Select **Checktune**.
4. Click  **Mass calibration/Checktune**.
5. Before the tune completes, click  **Stop Tune**.



5 Using Methods

The Method Editor Window	37
Set the Active Instrument Mode	38
Set Up and Run an Acquisition Method	39
Load the default method	39
Set Up a Scan Method	42
Run Methods	46
Review the Data using MassHunter Qualitative Analysis	49

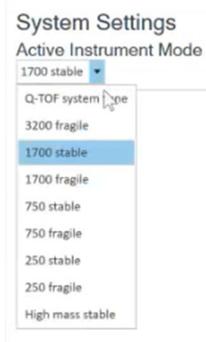
MassHunter Acquisition methods include the parameters for each component associated with your instrument.

The Method Editor Window

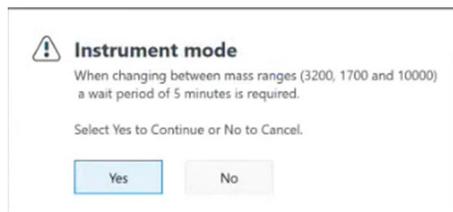
1. Launch the acquisition software: select **OpenLab Control Panel > Instruments** (bottom-left corner) **> your instrument > Launch**. Alternatively, if available double-click the desktop shortcut.
2. In the Windows section, select **Method Editor**. The Method Editor window opens in the Main Window.

Set the Active Instrument Mode

1. Click **Tune > Autotune** in the left pane.
2. Click  **Tune control**.
3. From the **Active Instrument Mode** drop-down list, select **1700 Stable**.



4. Select **Yes**. Wait for the system to stabilize.



5. Click  **Release tune control** in the toolbar to release control of the Q-TOF instrument.

Set Up and Run an Acquisition Method

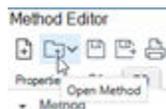


Working with the Default Method

Once an analysis has been created or opened, the default.m method is available to start from or apply a previously created method. The default method represents a good starting point for method development.

Load the default method

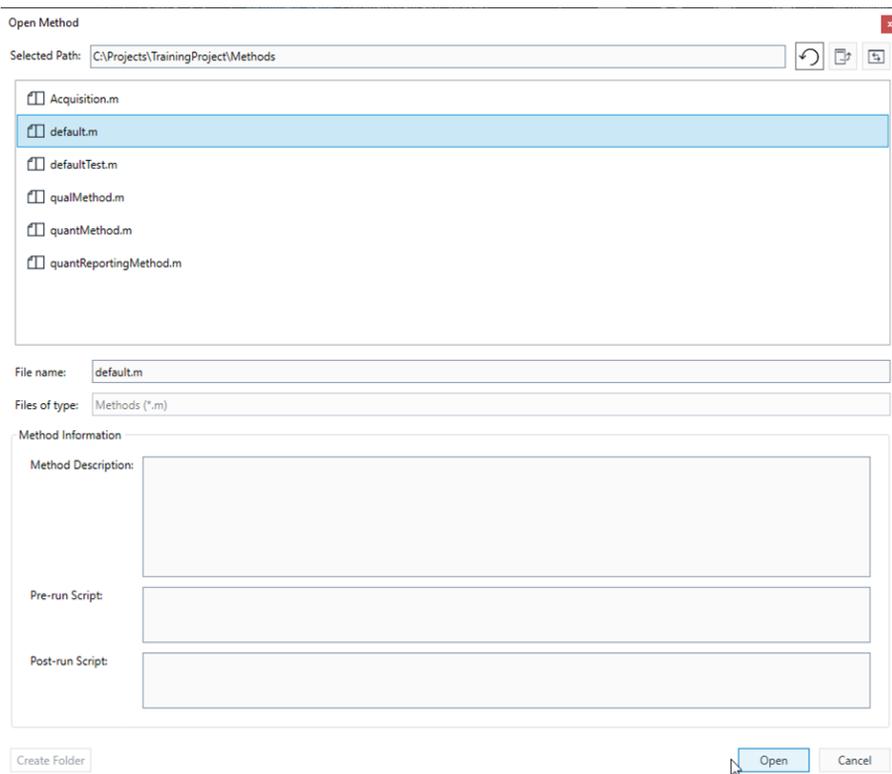
1. In the Method Editor window, click the **MS Q-TOF** tab.
2. Click **Open Method** to review the methods available.



Using Methods

Set Up and Run an Acquisition Method

3. Select **default.m** and click **Open**.



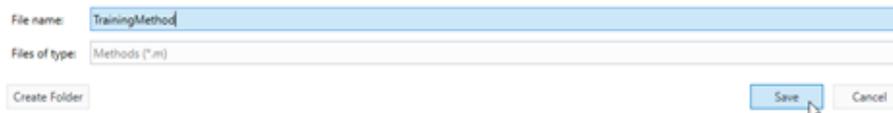
4. Under Method, review all the default Method subsections.
 - a. Acquisition – Set Q-TOF acquisition parameters.
 - b. Source – Set source parameters for the Q-TOF .
 - c. Chromatograms – Specify plots to display in the Chromatogram Plot window during the run.
5. Click **Save As Method**.



Using Methods

Set Up and Run an Acquisition Method

6. Enter a **File name**—for this example, *Training Method*—then click **Save**.



NOTE

After modifying or viewing a method using the drop-down list, you must apply the method to send the parameters to the instruments.

7. Click **Apply**.



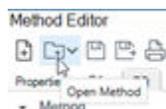
Set Up a Scan Method



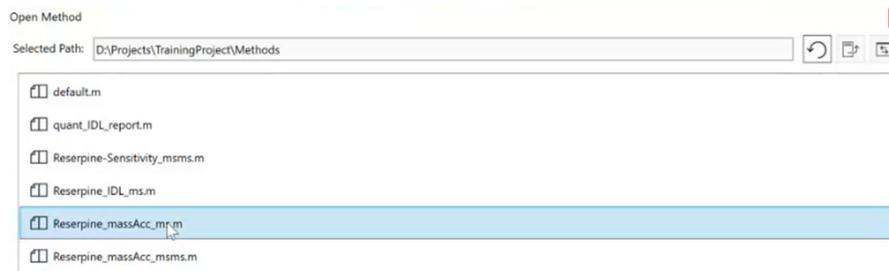
Load an existing method and Save As new method

In this exercise, you will use an existing method (the scan checkout method used during installation) to see if there is a signal for Reserpine at m/z 609 within the spectrum.

1. Click **Open Method** to review the methods available.



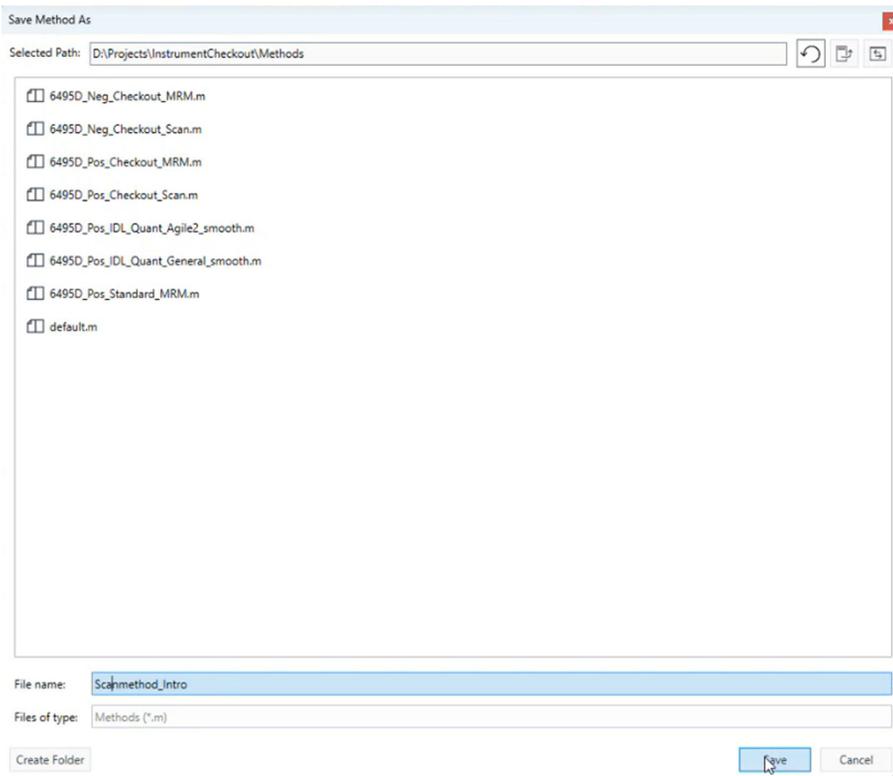
2. Select the checkout method used in installation (for example, Reserpine_massAcc_ms.m) and click **Open**. The method loads into the Method Editor.



Using Methods

Set Up a Scan Method

- Click **Save as Method**, enter *Scanmethod_Intro* for the file name, and click **Save**.



Using Methods

Set Up a Scan Method

- Select **Method > Acquisition**, then under Acquisition Parameters, review the acquisition settings, noting Start m/z and End m/z. Under TOF Acquisition Parameters, set the Rate to 3.

Acquisition Parameters

Polarity **Positive** Negative

MS Experiments

On	Start m/z	End m/z	Fragmentor (V)	Collision energy (V)	Skimmer (V)
1	100	1700	180	0	45

TOF Acquisition Parameters

MS-only Acquisition

Rate (spectra/s)

Time (ms/spectrum)

Transients/spectrum 5692

- Under MS Timetable, click **Add a row** at the end twice.

MS Timetable

Add row at the end

- Set the Start time (min) and Value as shown below.

Start time (min)	Type	Value
1	0 LC Stream Valve	To Waste
2	1 LC Stream Valve	To MS

- Click to enable **Post-run diverter position**.

Post-run diverter position To Waste

Using Methods

Set Up a Scan Method

- Click **Pump/Sampler/Column Comp settings** to review the settings programmed for the LC pump, noting the flow rate.

The screenshot displays the 'Binary Pump' configuration window. The 'Flow' section shows a rate of 0.400 mL/min. The 'Solvents' section is divided into two parts: the first part shows 100.0% Water V.03 for 0.20 minutes, and the second part shows 100.0% Acetonitrile V.03 for the rest of the run. The 'Pressure Limits' section has a minimum of 0.00 bar and a maximum of 600.00 bar. The 'Stoptime' is set to 4.35 minutes. The 'Advanced' section shows a 'Timetable (8/100 events)' table.

Time [min]	A [%]	B [%]	Flow [mL/min]	Max. Pressure Limit [bar]
0.00	50.00	50.00	0.400	600.00
0.20	50.00	50.00	0.400	---
0.85	10.00	90.00	0.400	---
2.85	10.00	90.00	0.400	---
2.86	50.00	50.00	0.400	---

- To send the current parameters shown in the Method Editor window to the LC and MS instruments, click **Apply**.

Run Methods



Run a Scan Method

In this exercise, you will acquire data using MassHunter Acquisition software and then use MassHunter Qualitative Analysis software to identify a precursor ion for Reserpine.

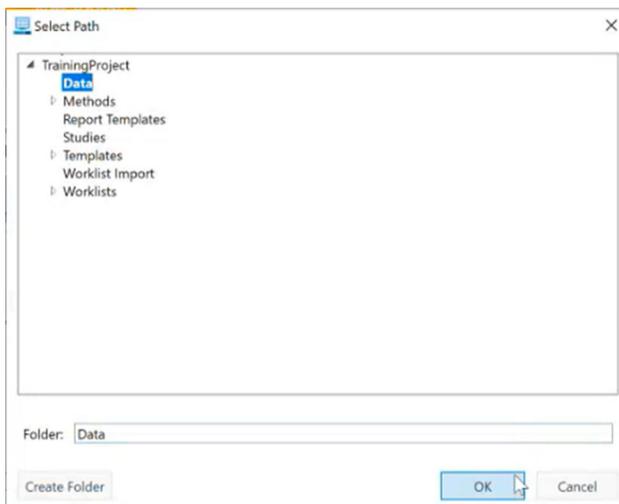
1. Place the Vial 4 100 pg/μl checkout sample, prepared during the system checkout, into a vial location of the sampler and note the location.
2. In the main window, click the **Sample Run** tab to display the Sample Run window.
3. In the Sample Run window, specify the following information:
 - a. Sample Name: Sample 1
Sample Position: Vial 4 (or applicable position)
 - b. Sample Injection Volume: Select **As Method** to use the volume specified in the method applied in the last step.
 - c. Data File Name: Introduction_Scan_001.d
(optional) Select **Auto Increment** to automatically increment the file name if that file exists.

NOTE

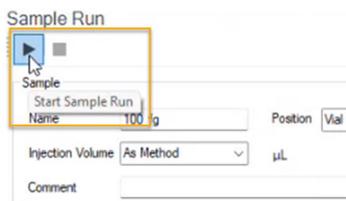
Using Methods

Run Methods

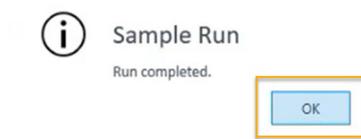
- d. Data File Path: Set to **D:\Projects\InstrumentCheckout\Data\Introduction**. Create a folder if necessary.



4. Click **Start Sample Run**.



5. Click **OK** when the run completes.



**Monitor the Status Windows**

As data is being acquired, use the instrument status monitors and online signal displays available in the Instrument Status and Real-time Plot Panes to observe changes in modules.

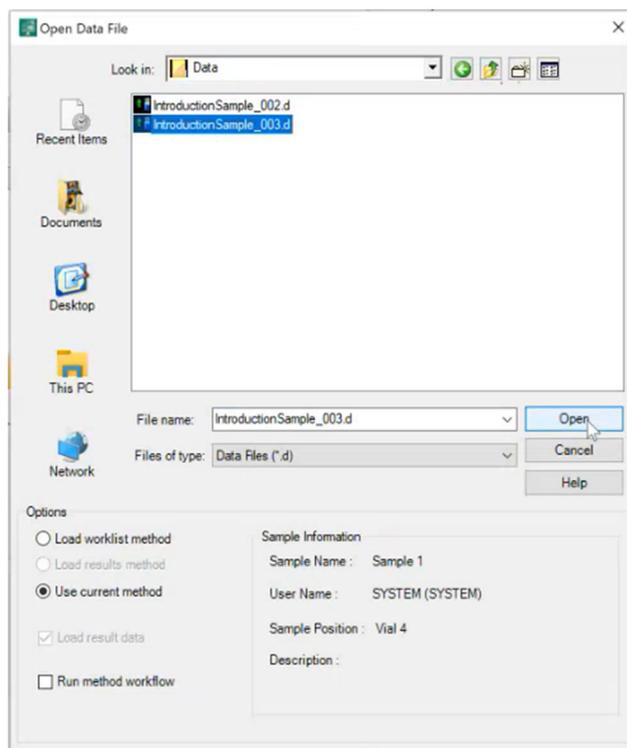
1. View the Chromatogram Plot and note the retention time for Reserpine.
2. Observe the Spectrum window while the samples runs. Discuss with your Agilent-certified service professional the changes observed over time.

Review the Data using MassHunter Qualitative Analysis

1. From the desktop, double-click the MassHunter Qualitative Analysis icon.



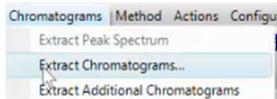
2. In the Open Data File window, browse to the data file directory created earlier (Introduction), select the data file to review, and click **Open**.



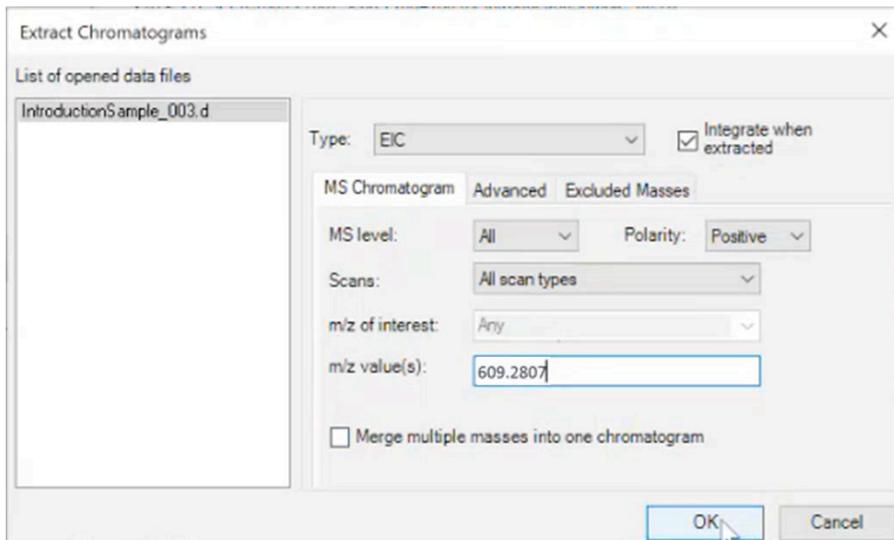
Using Methods

Review the Data using MassHunter Qualitative Analysis

3. In the main window, click **Chromatograms > Extract Chromatograms**.



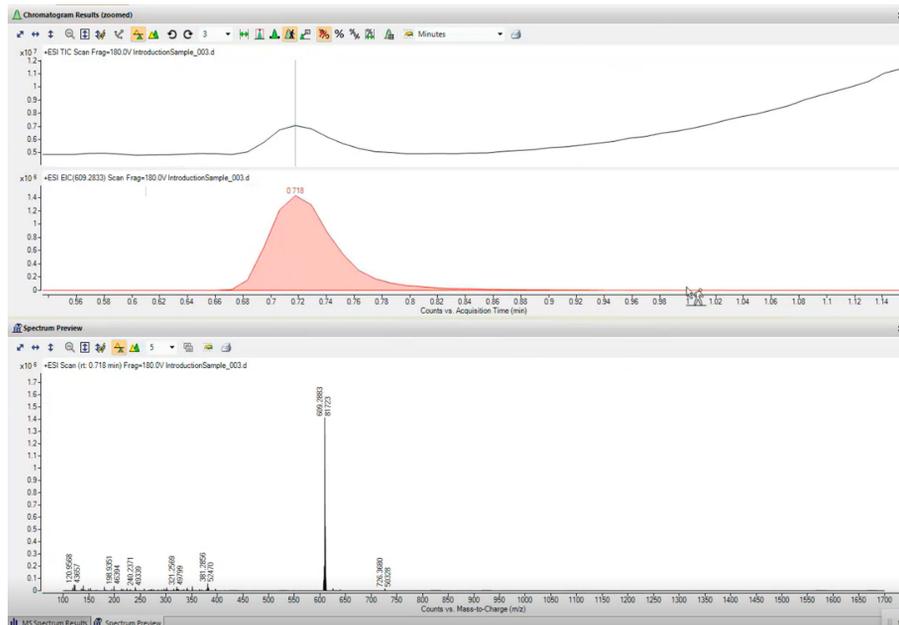
4. In the Extract Chromatograms dialog box, click **Type:** and select **EIC**.
5. Enter the *m/z* value: 609.2807, then click **OK**.



Using Methods

Review the Data using MassHunter Qualitative Analysis

6. Review the results.



Review the Results

1. What is the retention time for Reserpine?

2. What is the m/z observed for Reserpine in the mass spectrum?



6 Run a Worklist

Create and edit a worklist	53
Add Multiple Samples	54
Change Result Window Layouts	56
Use layout icons on the toolbar	56
Change the panes in the Compound Information window for Cal-L4	57
Save the default layout without the calibration curve	58
Load the newly created layout	59
Recreate (do not restore) the default layout	61

Run a Worklist

Create and edit a worklist

Use the Worklist window to create a list of samples to run. A worklist can be created using either the Study Manager Program or the Worklist window. Use this procedure to inject multiple samples by creating a new worklist.

Create and edit a worklist

1. In MassHunter Acquisition, click **Worklist** to show the Worklist window.



2. In the Worklist window, click **Open Worklist**.
3. Select the G6530_DAJ.S.w worklist and click **Open**.
4. In the Worklist window, click  **Add Multiple Samples**. The Add Multiple Samples dialog opens.

NOTE

Samples can also be added one-by-one (for example, if you only need to run a few samples or several replicates of the same sample).

Add Multiple Samples

1. On the **Sample Position** tab, specify the sample vial locations (make sure the specific sample tray type has been configured by right clicking the autosampler device image). Click **OK**.

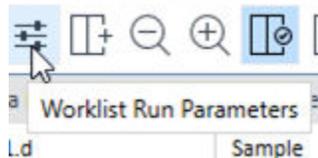
The screenshot shows the 'Add Multiple Samples' dialog box with the 'Sample Position' tab selected. The dialog is divided into several sections:

- Sample Information:** Includes a 'Sample Position' tab and a 'Current Configuration' checkbox.
- Select Well-plate or Vial Tray:** A tree view showing 'Well-plate Tray' expanded to show 'Plate 1' through 'Plate 8'.
- Plate/Tray Type:** A dropdown menu set to '6 Vials Generic Plate'.
- Selection Origin:** Radio buttons for 'Top left' (selected), 'Top right', 'Bottom left', and 'Bottom right'.
- Block Increment:** Radio buttons for 'Row major' (selected), 'Column major', and 'Serpentine'.
- Number of samples:** A text box containing '0'.
- Number of replicates:** A text box containing '1'.
- Well-plate/Tray:** A 2x3 grid of vial positions. The grid is labeled with columns 1, 2, 3 and rows A, B. The entire grid is highlighted with a red border.
- Buttons:** 'OK' and 'Cancel' buttons at the bottom right.

Run a Worklist

Add Multiple Samples

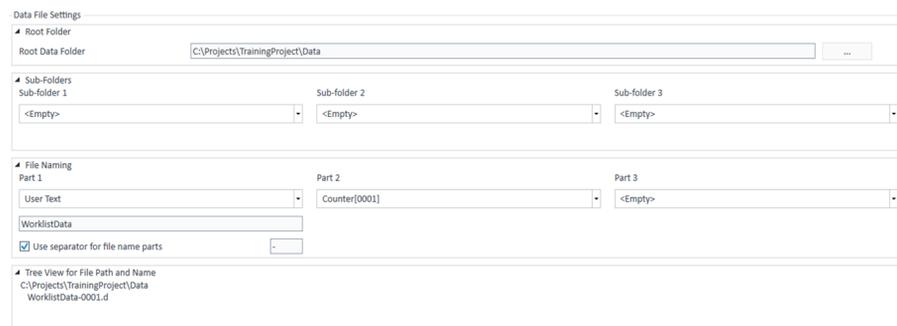
- To set up the worklist run, click **Worklist Run Parameters**.



- On the **Run Parameters** tab, type the paths for the method.



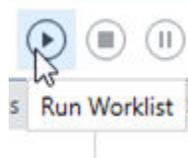
- In the **Data File Settings** section, expand all options and enter or select the folders for the data files. Select the file naming options. Click **OK**.



NOTE

For more information on the **Intelligent Reflex** tab, see the online Help.

- Optional. On the **Additional Parameters** tab, enter a comment, and click **OK**.
- To start the worklist, click **Run Worklist**.

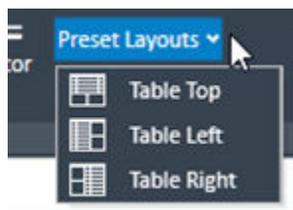


Change Result Window Layouts

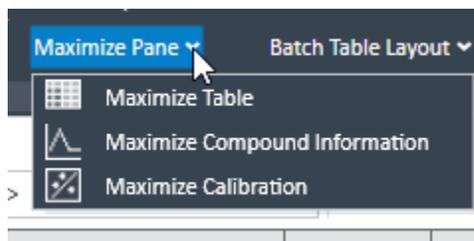
This task shows you how to customize your layout and how to recreate the default layout.

Use layout icons on the toolbar

1. Use layout options on the View Tab to position the Batch Table, Compound Information, and Calibration Curve windows.
 - a. On the **View** tab, select **Preset Layouts > Table Left**.
 - b. On the **View** tab, select **Preset Layouts > Table Right**.
 - c. On the **View** tab, select **Preset Layouts > Table Top**.



2. Use layout icons on the View Tab to maximize each individual window:
 - a. On the **View** tab, select **Maximize Pane > Maximize Table**.
 - b. On the **View** tab, select **Maximize Pane > Maximize Compound Information**.
 - c. On the **View** tab, select **Maximize Pane > Maximize Calibration**.



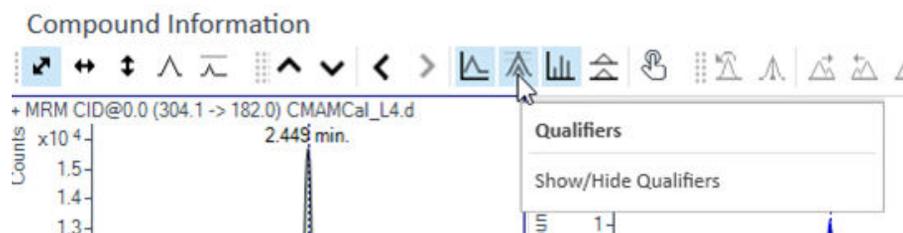
Run a Worklist

Change Result Window Layouts

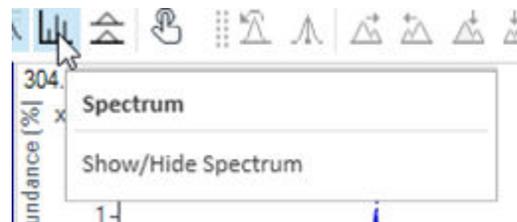
3. On the **View** tab, click **Restore Default Layout**.

Change the panes in the Compound Information window for Cal-L4

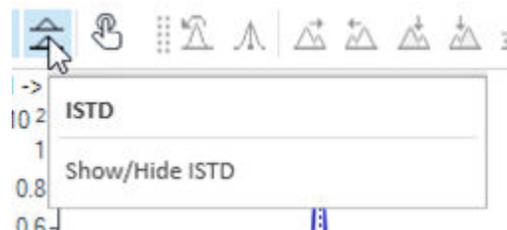
1. In the Batch Table, select the **Cal-L4** row.
2. On the Compound Information toolbar, click **Show/Hide Qualifiers**.



3. Click **Show/Hide Spectrum**.



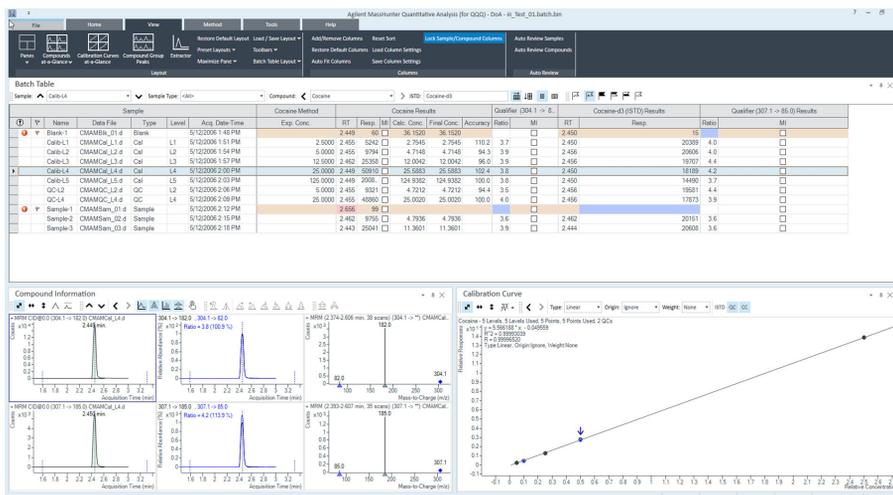
4. Click **Show/Hide ISTD**.



Run a Worklist

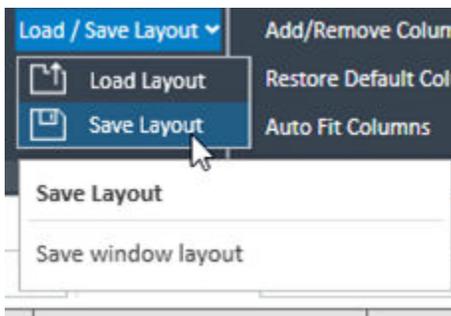
Change Result Window Layouts

- The layout and results look like those in the following figure.



Save the default layout without the calibration curve

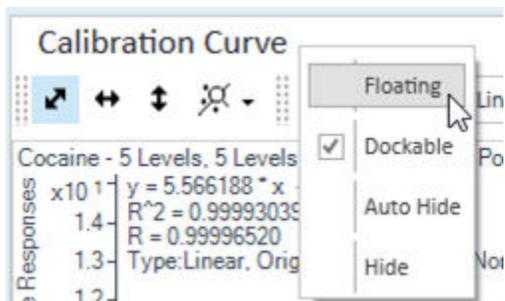
- Close the Calibration Curve window.
- On the **View** tab, select **Load/Save Layout > Save Layout**.



- In the Save Layout File dialog, name the layout file *Batch Table plus Compound Information* and click **Save**.

Load the newly created layout

1. On the **View** tab, click **Restore Default Layout**.
2. On the **View** tab, select **Load/Save Layout > Load Layout**.
3. In the Load Layout dialog, select **Batch Table plus Compound Information** and click **Open**.
4. On the **View** tab, click **Restore Default Layout**.
5. Right-click inside the title bar of the Calibration Curve window, and select the **Floating** check box.



6. Right-click the title bar of the Compound Information window, and select **Floating**.

Run a Worklist

Change Result Window Layouts

10. Right-click inside the title bar of the Calibration Curve window and uncheck the **Floating** check box.
11. Move the Compound Information window so that the layout corresponds to the one pictured at the start of the task.

Recreate (do not restore) the default layout

1. Maximize the program main view.
 - Anchor the Calibration Curve window first, and then the Compound Information window, to recreate the default layout.
 - If after anchoring the two windows, the calibration curve is on the left side, right-click the title bar of the Calibration Curve window and drag it to the right. A gray rectangle shows where this window will be placed within the main view.
 - Drag the calibration curve to the bottom-right corner of the main view.
2. On the **View** tab, click **Restore Default Layout**.



7

Maintenance

Register on Agilent SubscribeNet (New Account Registration)	63
Perform Back Up and Platform Best Practices	64

Register on Agilent SubscribeNet (New Account Registration)

NOTE

If you already have a SubscribeNet account set up for previously purchased Agilent products, it is not necessary to set up additional accounts.

1. Using a web browser, navigate to <https://agilent.subscribenet.com/>.

Agilent Technologies Agilent SubscribeNet

Electronic Software and License Delivery

Please login. Your Login ID is your Email address.

Login ID

Password

Remember my password until I logout

Login

If you have forgotten your login ID, password, or are not sure whether you have an account use our [Password Finder](#).

SubscribeNet new account registration.

Customers who have an authorization code from their Agilent product purchase may [CLICK HERE](#) to register and create a new SubscribeNet Account and Login ID.

[Privacy Statement](#) | [Terms of Use](#) | [Agilent Home](#) | © Agilent 2000-2023

2. Click **CLICK HERE** to register.
3. Enter the following required information to create the account, along with the authorization code received from the product purchase.
 - a. Authorization Code
 - b. Email
 - c. Company
 - d. Department

Maintenance

Perform Back Up and Platform Best Practices

- e. First Name
 - f. Last Name
 - g. Phone
 - h. Address
 - i. City
 - j. State/Province
 - k. Country
4. Click **Submit**.
You'll receive an email to activate your account.

Perform Back Up and Platform Best Practices

- Safely store the software media provided for the system.
- Set up data/computer image backups regularly.
 - [Microsoft Back up and Restore Options](https://t.ly/r2995) (https://t.ly/r2995)
- Disable power management options and automatic utilities.
 - Set power options to **Put the computer to sleep = Never**.
 - Set Windows Update to **Check for updates but let me choose whether to download and install them**.
- Turn on Windows Firewall.
 - Select **Notify me when Windows Firewall block a new program**.

Maintenance

Perform Back Up and Platform Best Practices



Locate the tune solution bottle and properly store tune solution

Locate the tune solution bottle and confirm that it is safely stored in the appropriate temperature conditions.



Perform daily cleaning of Ionization Source and Spray Chamber

Perform this maintenance daily or at the end of each shift or anytime you suspect carryover contamination from one sample or analysis to another.

After determining the Source type, find the instructions for cleaning in the user guide for the source in use.



Review routine procedures in the user guide

Perform maintenance daily or at the end of each shift or anytime you suspect carryover contamination from one sample or analysis to another.



Remove, clean, and replace the capillary

Review the steps in the user guide.



Add a new user defined EMF counter

Using the online help, set up a user defined EMF counter by entering a new threshold value for a selected EMF item.

www.agilent.com

© Agilent Technologies, Inc.

G1960-90153

January 2026

