



6545XT LC/Q-TOF System **Introduction Workbook**

with OpenLab CDS Software



Notices

Document Identification

G1960-90158

May 2026

Copyright

© Agilent Technologies, Inc. 2026

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This workbook is valid for OpenLab CDS Acquisition software 3.0, until superseded.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	8
	Additional resources	8
	Where to find more information	10
2	Hardware	11
	Overview	12
	Front view	13
	Side view	14
	Back view	15
	Basic Components	16
	Ionization source	16
	Calibrant delivery system	18
	Nebulizer	19
	Rough pump	20
	Waste bottle from LC pump or mass spectrometer	21
3	Software Basics	22
	Software Start-Up	24
	User Interface and General Navigation	25
	Close Connection	26
	Creating and Configuring Projects	27
	Instruments	29
	Create an instrument	30
	Launch an instrument	31
	Create an instrument shortcut	32
4	Tuning	33
	Mass Calibration	35
	Checktune	37
	Generating a Detailed Tune Report	39

5	Using Methods	41
	The Method Editor Window	43
	Set the Active Instrument Mode	44
	Set Up and Run an Acquisition Method	45
	Load the default method	45
	Load an Acquisition Method	47
	Run Methods	51
	Review the Data using MassHunter Qualitative Analysis	54
6	Setting Up and Running a Sequence	57
7	Maintenance	61
	Register on Agilent SubscribeNet (New Account Registration)	62
	Perform Back Up and Platform Best Practices	64



1

Introduction

About this Training Workbook	6
How to use this Workbook	7
Before You Begin	8
Additional resources	8
Where to find more information	10

About this Training Workbook

This workbook provides instructions for the Agilent 6545XT LC/Q-TOF system running OpenLab Acquisition 3.0 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 6545XT AdvanceBio 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 6545XT AdvanceBio Q-TOF LC/MS System User Guide* by scanning the code or navigating to <https://aglt.co/LCMSUserDocs>.
- Use the *Agilent 6545XT AdvanceBio 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>.

Agilent Community

Introduction

Before You Begin



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/>

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	12
Front view	13
Side view	14
Back view	15
Basic Components	16
Ionization source	16
Calibrant delivery system	18
Nebulizer	19
Rough pump	20
Waste bottle from LC pump or mass spectrometer	21

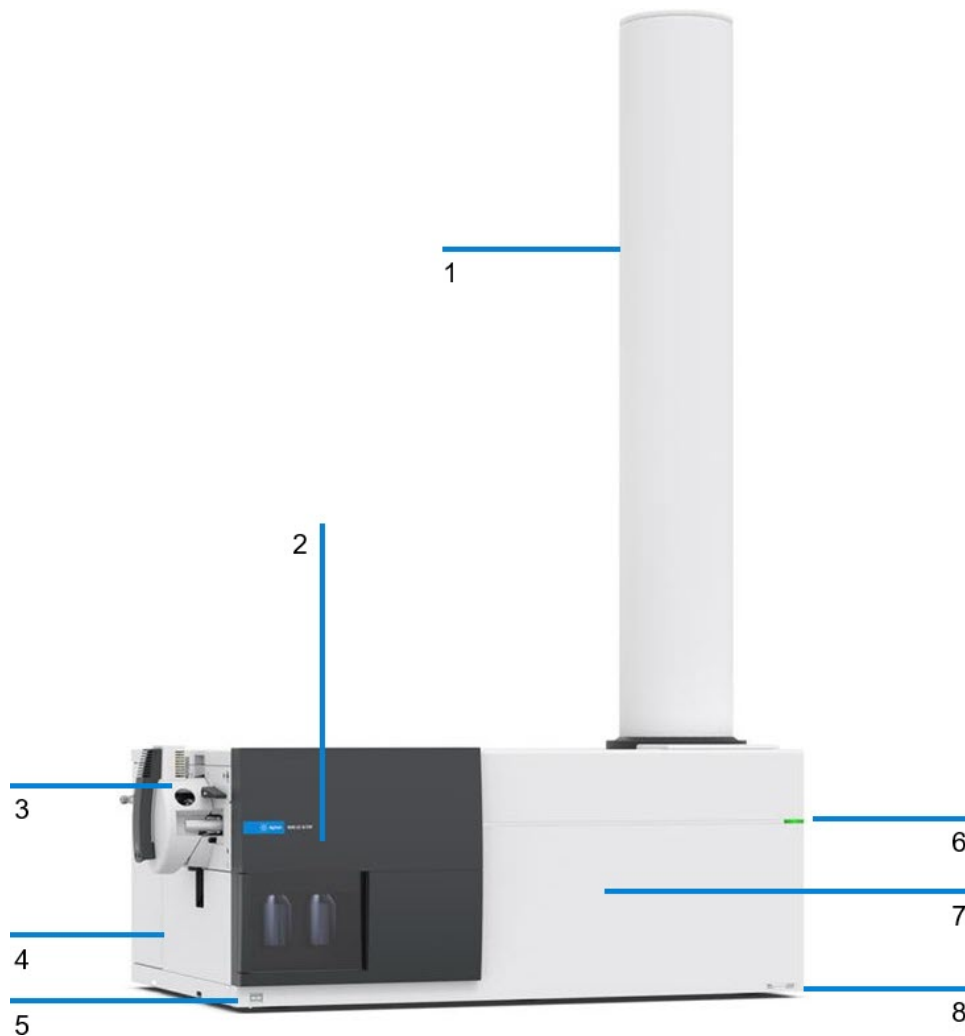
Overview

In this section, you will identify basic hardware components and their locations for the Agilent 6545XT AdvanceBio LC/Q-TOF system.

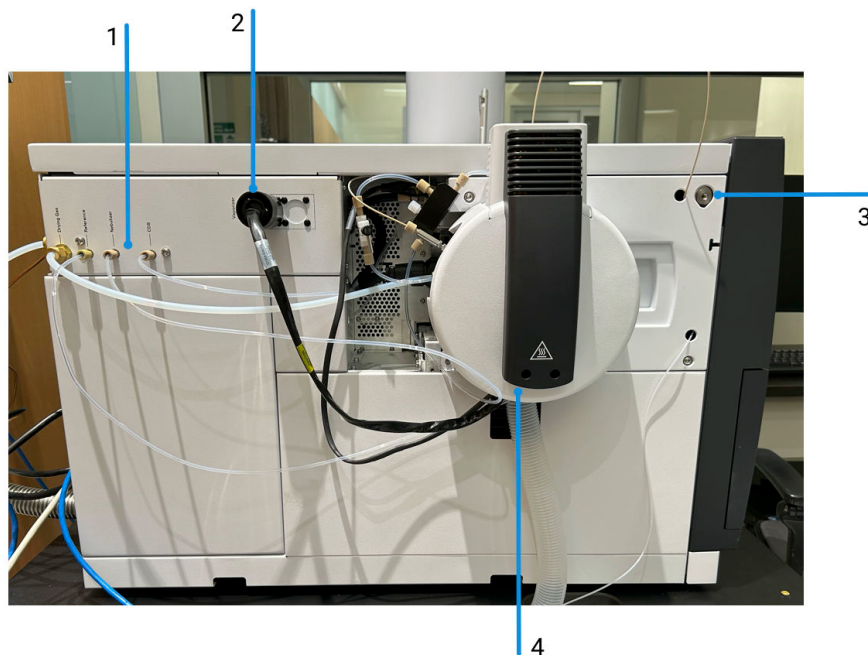


Fill in the Blank

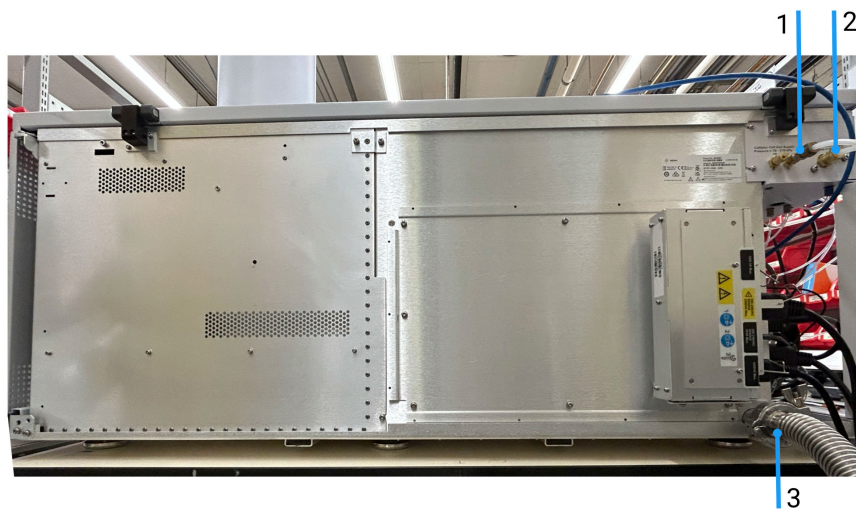
Fill in the Blank: Work with your Agilent Service Engineer and/or use the *Agilent 6545XT AdvanceBio Q-TOF LC/MS System User Guide* to label the flagged components below for your installed instrument(s).

Front view

Side view



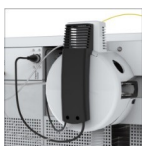
Back view



Basic Components

Ionization source

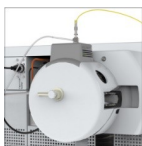
Agilent liquid chromatography/mass spectrometry (LC/MS) ion sources enable analysis of a wide range of samples quickly and accurately. These easily interchangeable ion sources for Agilent LC/MS systems include the:



- Agilent Jet Stream (Dual AJS ESI)



- Electrospray Ionization (ESI)



- Multimode Ionization source (MMI)



- Agilent nanospray ESI

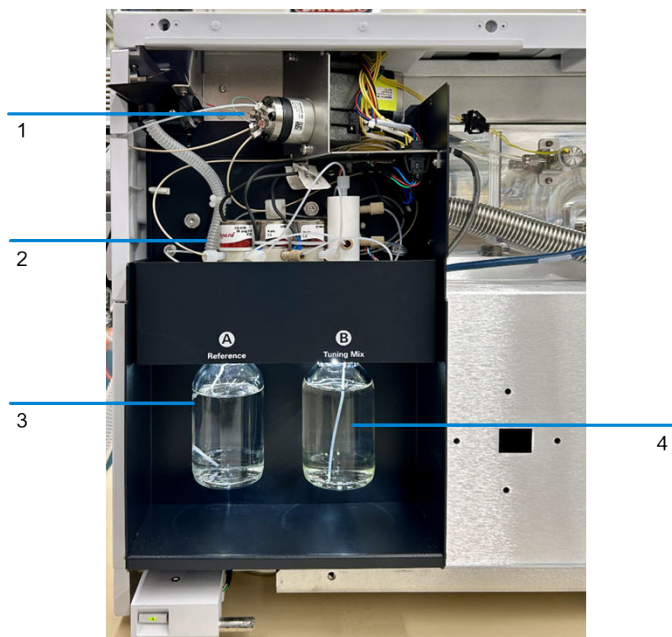
**Hardware Introduction**

1. List the ionization source in use:

2. 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.



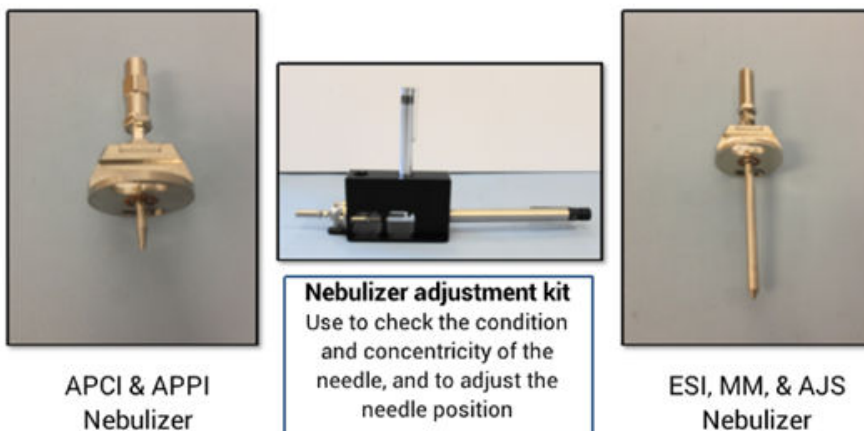
Hardware Introduction

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+



For more information on your rough pump, visit <https://www.agilent.com/en/product/vacuum-technologies>.

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	24
User Interface and General Navigation	25
Close Connection	26
Creating and Configuring Projects	27
Instruments	29
Create an instrument	30
Launch an instrument	31
Create an instrument shortcut	32

The OpenLab Control Panel is the administrative and management center for the OpenLab 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

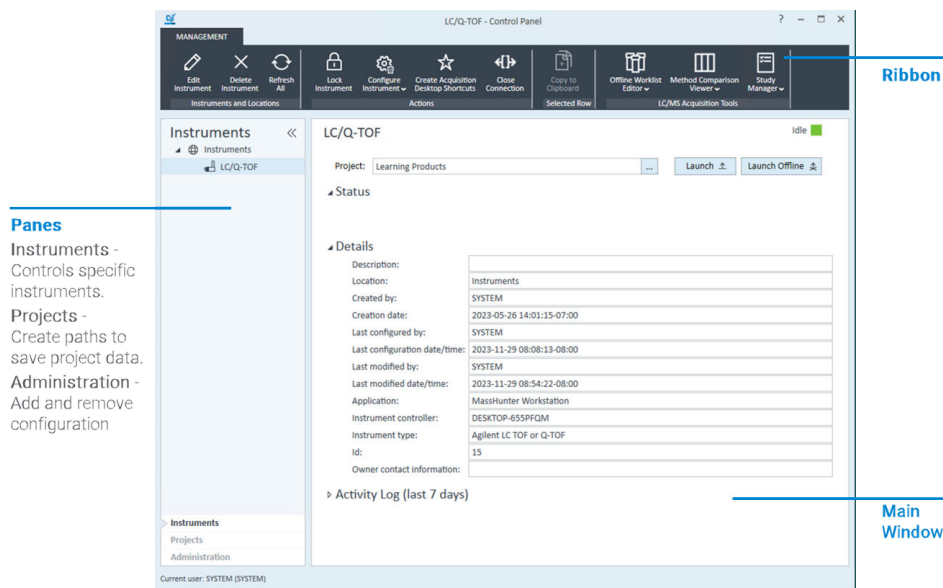


- From the desktop, double-click the **Control Panel** icon.



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

User Interface and General Navigation



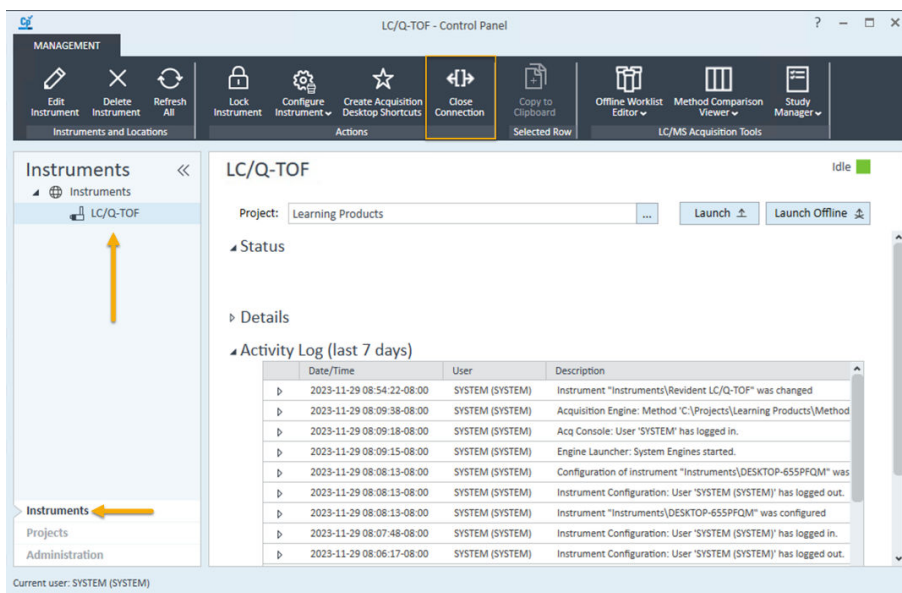
Adjusting the Instruments pane:

- 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.

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

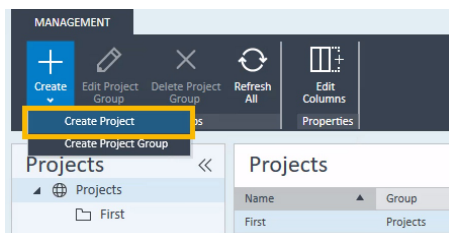


Creating and Configuring Projects



Creating and Configuring Projects

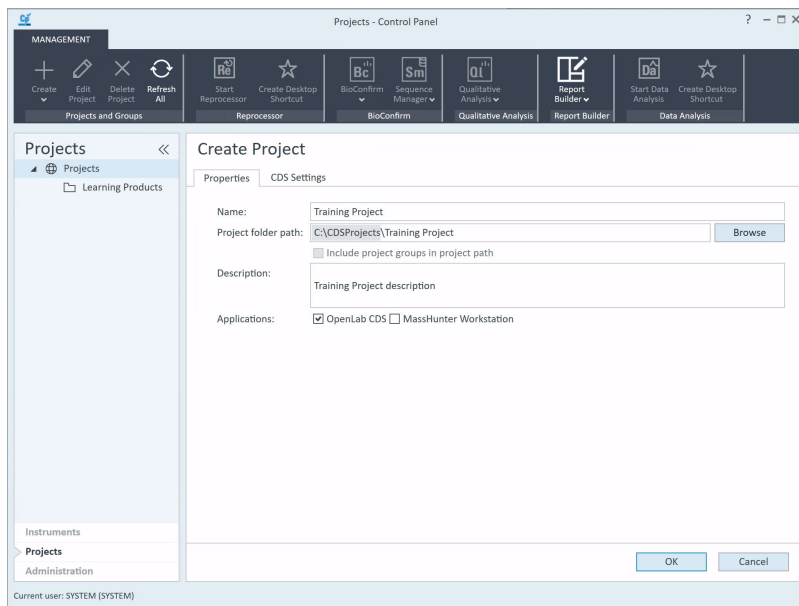
1. Click  **Projects** and select **Projects**.
2. On the ribbon, click **Create Project**.



3. In the **Name** text box, type *Training Project*.
4. In the Project folder path text box, leave the default folder path.
5. (optional) In the Description text box, type a description of the project, for this example *Training Project Description*.
6. If necessary, select the **OpenLab CDS** checkbox.
7. Click the **CDS Settings** tab and review the available options. Do not change the defaults.
8. Click **OK**.

Software Basics

Creating and Configuring Projects



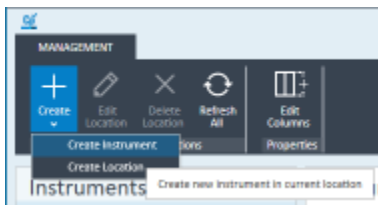
Instruments

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**.

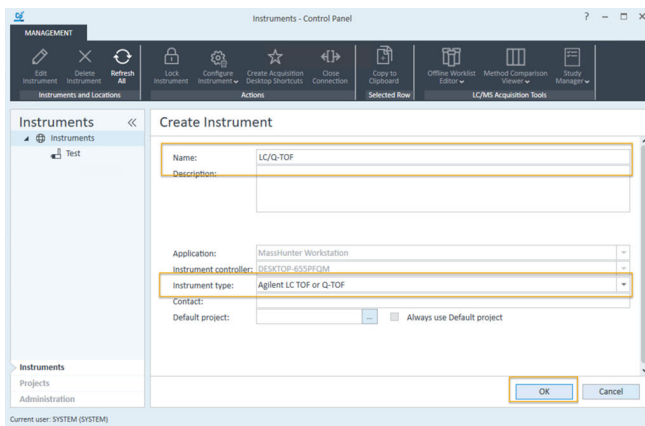


3. Enter the required data in the Create Instrument pane.
 - a. Name: 6545XT AdvanceBio LC/Q-TOF
 - b. Instrument Type: **Agilent LC TOF or Q-TOF**

NOTE

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

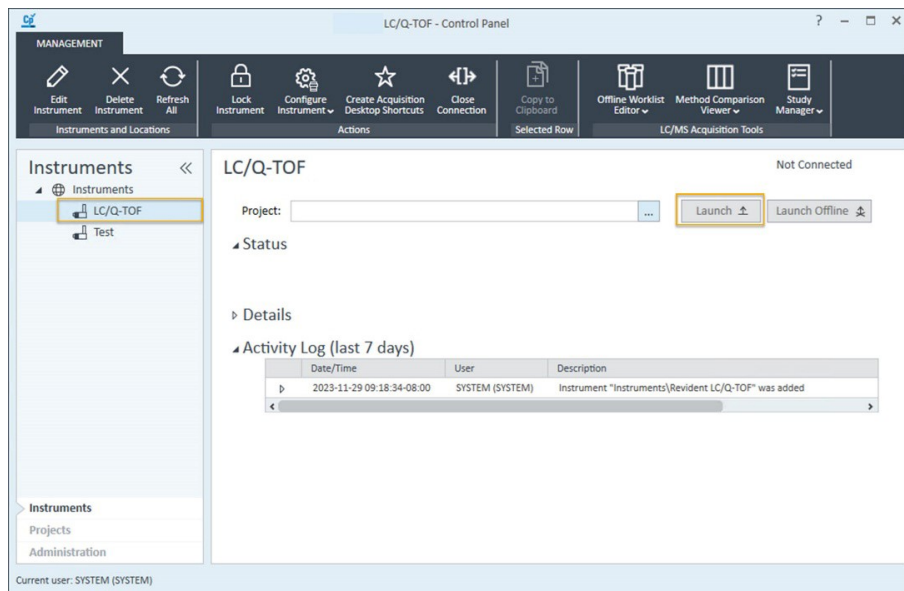
4. Click **OK**.
5. Click ... to select the **TrainingProject** project from the Select Project dialog.
6. Click **OK**. The instrument is displayed in the navigation pane.



Launch an instrument

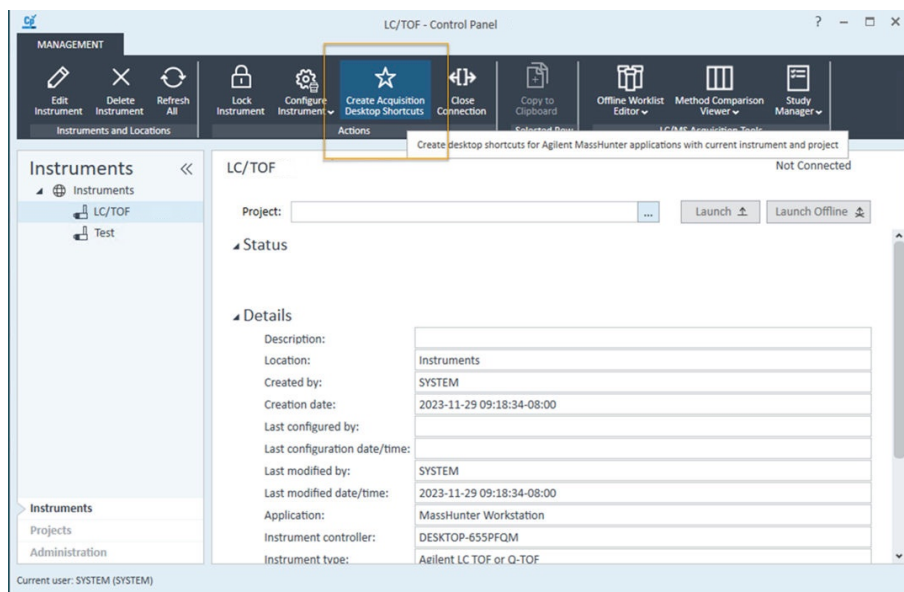
To begin acquisition, launch the instrument from the instrument table or the instrument details page, or launch an instrument directly from your desktop shortcut.

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



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

Mass Calibration	35
Checktune	37
Generating a Detailed Tune Report	39

In this chapter, you will run a mass calibration and a checktune, and generate a tune report.

Mass Calibration

Mass Calibration performs a TOF mass calibration on the active instrument mode by infusing a solution with known masses into the source and measuring the actual flight times for ions of the masses.

For optimum performance, run a mass calibration before analysis each day, or after changing high-voltage manual tune parameters or adjusting the mass range. During normal operation, mass calibration takes approximately 5 minutes. To ensure the mass accuracy of your data, a mass calibration is required when selecting a different active instrument mode.

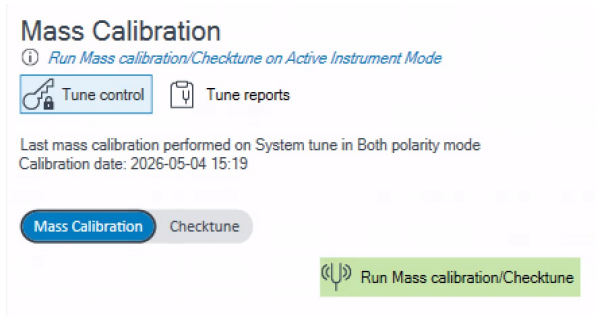
NOTE

Only the active instrument mode can be calibrated.

A mass calibration can only be run on instrument modes that have first been configured and enabled in the Configure Tune panel. System Tune is always enabled by default.

Follow these steps to perform a mass calibration:

1. In the Method Editor window, click the **MS TOF/Q-TOF** tab.
2. Click **Tune > Mass Calibration** in the left pane.
3. Click  **Tune control**.
4. Select **Mass Calibration**.



Tuning

Mass Calibration

5. Click  **Mass calibration/Checktune**. When the tune completes, review the report.
6. Click  **Tune control** to release control of the LC/Q-TOF instrument.

Checktune

Checktune is used to determine whether the instrument is correctly tuned by applying the current tune parameters to the source, quadrupole, and detector to confirm that the signal intensity and resolution is within the expected range. A checktune performs a TOF mass calibration, checks the Quad, and checks the detector for the currently selected active instrument mode. A checktune is automatically performed when running an autotune.

During normal operation, a checktune will take approximately 20 minutes.

A checktune can be run with the following ion sources:

- ESI
- Dual AJS ESI

NOTE

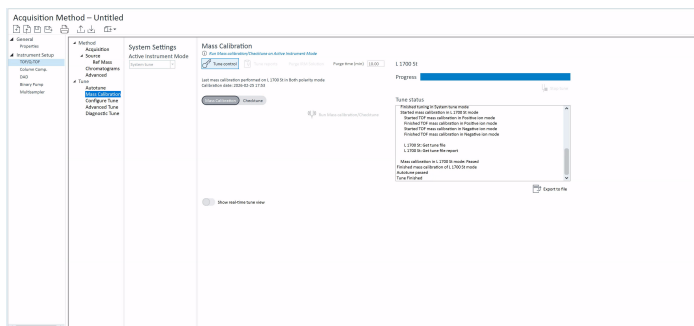
A checktune is only run on the currently selected instrument mode in the Active Instrument Mode drop-down.

A Checktune can only be run on instrument modes that have first been configured and enabled in the Configure Tune panel. System Tune is always enabled by default.

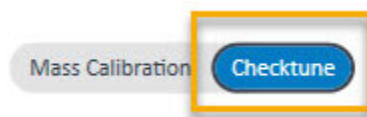
Follow these steps to perform a checktune:

1. In the Method Editor window, click the **MS TOF/Q-TOF** tab.
2. Click **Tune > Mass Calibration** in the left pane.

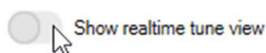
3. Click  **Tune control** **Tune control**.



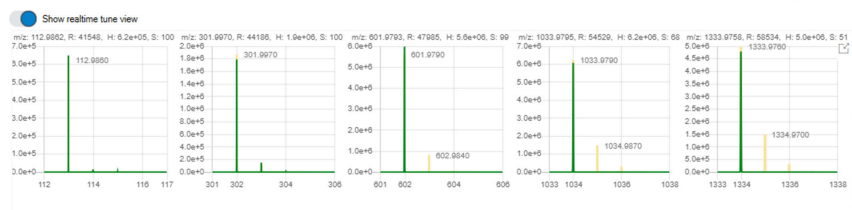
4. Select **Checktune**.



5. Activate **Show realtime tune view** to view the real time data acquisition.



6. Click  **Run Mass calibration/Checktune**. View the progress of the tune in the Real Time tune results.

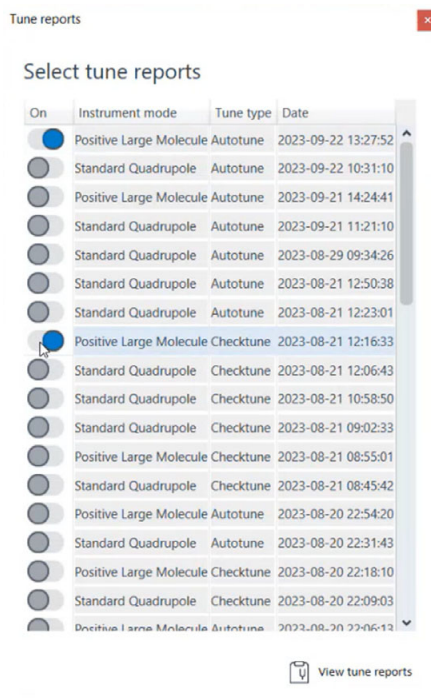


7. When the tune completes, review the report.
8. Click  **Tune control** **Tune control** to release control of the LC/Q-TOF instrument.

Generating a Detailed Tune Report

Generate a detailed tune report after running an Autotune, Checktune, or Mass Calibration.

1. In the Method Editor window, click the **MS TOF/Q-TOF** tab.
2. Click **Tune > Autotune** in the left pane.
3. Click  **Tune control**.
4. Click  **Tune Reports**. If this button is not available, run an autotune, checktune, or mass calibration first.
5. Use the toggles to select reports for viewing.



Tuning

Generating a Detailed Tune Report

6. Click  **Tune Reports** to generate PDFs for viewing in a PDF viewer.
7. Click  **Tune control** to release control of the Q-TOF instrument.



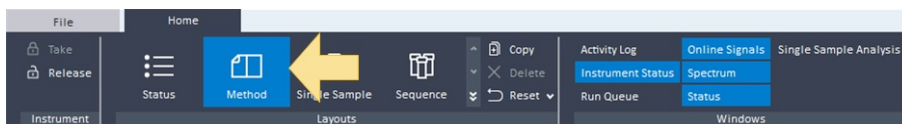
5 Using Methods

The Method Editor Window	43
Set the Active Instrument Mode	44
Set Up and Run an Acquisition Method	45
Load the default method	45
Load an Acquisition Method	47
Run Methods	51
Review the Data using MassHunter Qualitative Analysis	54

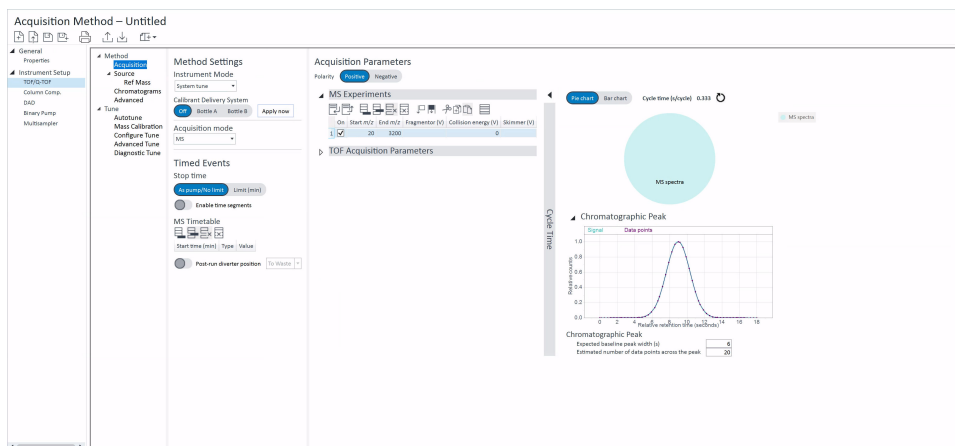
OpenLab Acquisition methods include the parameters for each component associated with your instrument.

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 OpenLab Acquisition, click **Method**.

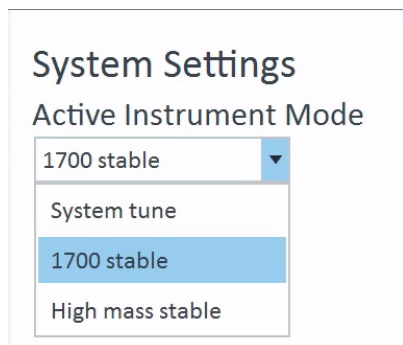


The Method Editor opens in the Main Window.

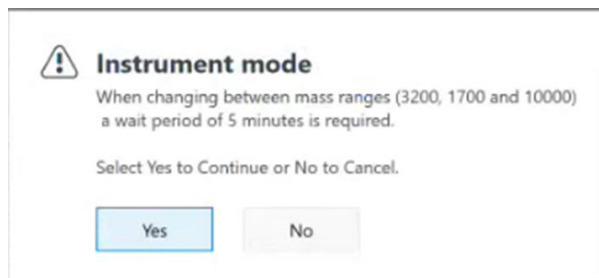


Set the Active Instrument Mode

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

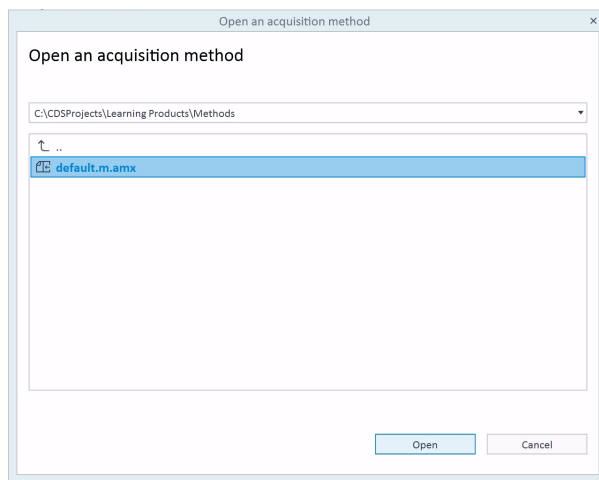


Working with the Default Method

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  **Open Method** to review available methods.
2. Select **default.m** and click **Open**.

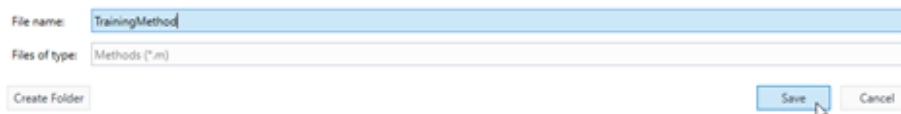


3. Under Method, review all the default Method subsections.
 - a. Acquisition – Set Q-TOF acquisition parameters.
 - b. Source – Set source parameters for the Q-TOF.
 - c. Chromatograms – Specify plots to display in the Chromatogram Plot window during the run.

Using Methods

Set Up and Run an Acquisition Method

- d. Advanced – Set advanced parameters such as Data Storage Thresholds, Data Storage Mode, and Experiment Defaults.
4. Click  **Save As Method**.
5. Enter a **File name**—for this example, *Training Method*—then click **Save**.



File name: TrainingMethod

Files of type: Methods (*.m)

Create Folder Save Cancel

NOTE

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

6. Click  **Apply**.

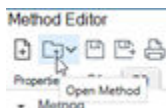
Load an Acquisition Method



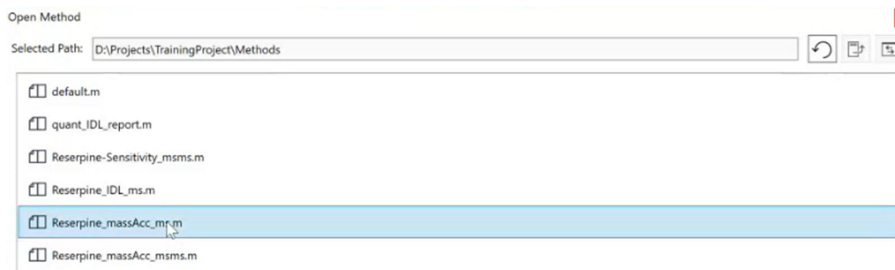
Load an existing method and Save As new method

In this exercise, you will use an existing method (the mass accuracy 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.



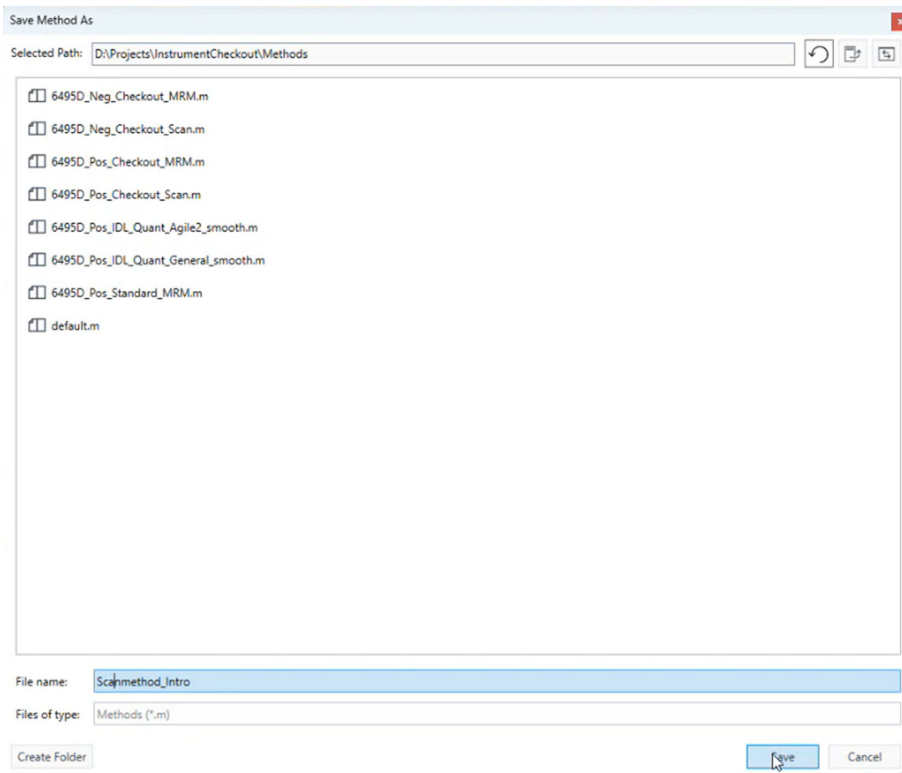
2. Select the checkout method used in installation (for example, Reserpine_massAcc_ms.m) and click **Open**. The method loads into the Method Editor.



Using Methods

Load an Acquisition Method

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



Using Methods

Load an Acquisition Method

4. 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

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

MS Timetable

6. 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

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

☒ Post-run diverter position

Using Methods

Load an Acquisition Method

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

Properties DA Multisampler Multisampler Pretreatment **Binary Pump** Column Comp MS Q-TOF

Flow: 0.400 mL/min

Solvents:

1	2
100.0 % Water V.03	100.0 % Water V.03
100.0 % Acetonitrile V.03	100.0 % Acetonitrile V.03

Pressure Limits: Min: 0.00 bar, Max: 600.00 bar

Stoptime: 4.35 min

Advanced: Timetable (8/100 events)

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	---

Buttons: Add, Remove, Clear All, Clear Empty, Cut, Copy, Paste, Shift Times

ISFT

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

Run Methods



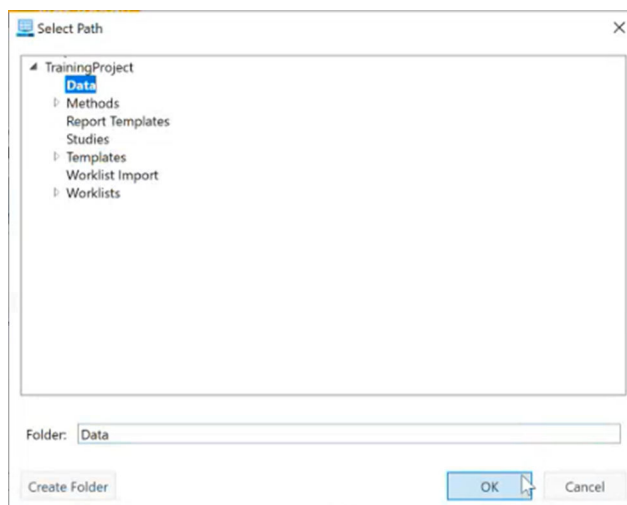
Run an acquisition method

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

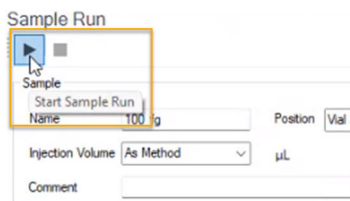
1. Place the Vial 4 100 pg/μl checkout sample, prepared during the system checkout, into a vial location of the sampler and note the location.
2. In the main window, click the **Sample Run** tab to display the Sample Run window.
3. In the Sample Run window, specify the following information:
 - a. Sample 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_001.sirslr
(optional) Select **Auto Increment** to automatically increment the file name if that file exists.

NOTE

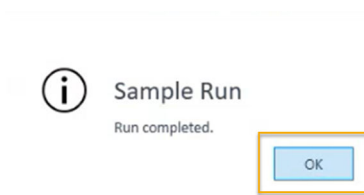
- d. Data File Path: Set to **D:\Projects\InstrumentCheckout\Data\Introduction**. Create a folder if necessary.



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.

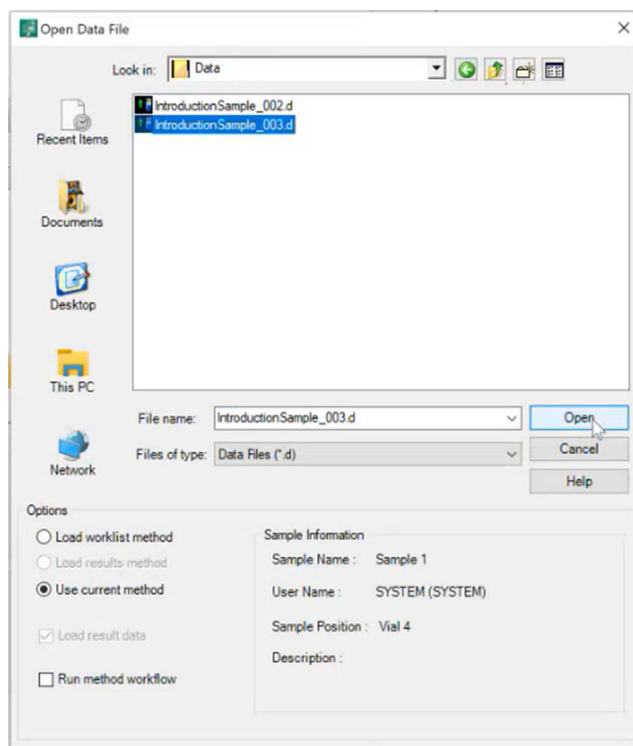
1. View the Chromatogram Plot and note the retention time for Reserpine.
2. Observe the Spectrum window while the samples runs. Discuss with your Agilent-certified service professional the changes observed over time.

Review the Data using MassHunter Qualitative Analysis

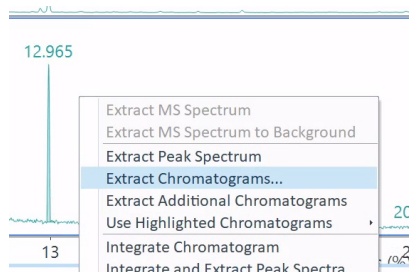
1. From the desktop, double-click the MassHunter Qualitative Analysis icon.



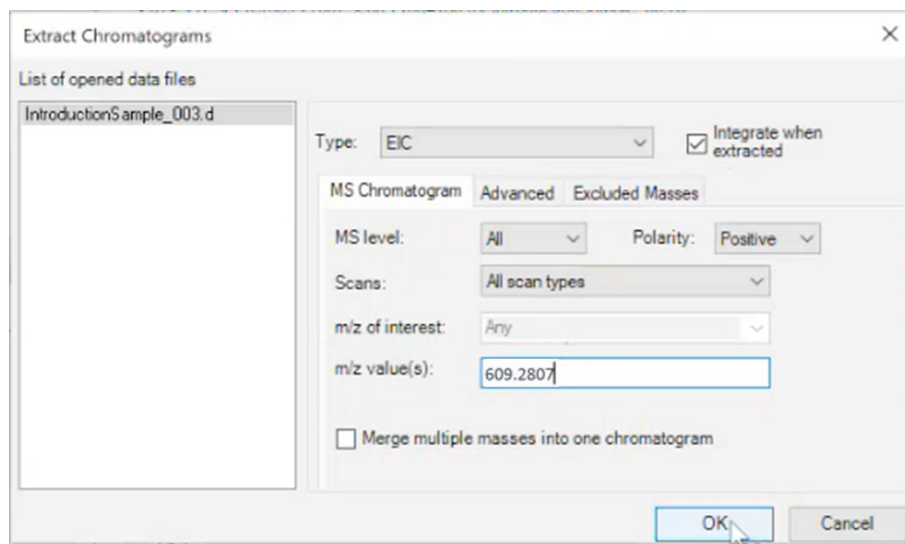
2. In the Open Data File window, browse to the data file directory created earlier (Introduction), select the data file to review, and click **Open**.



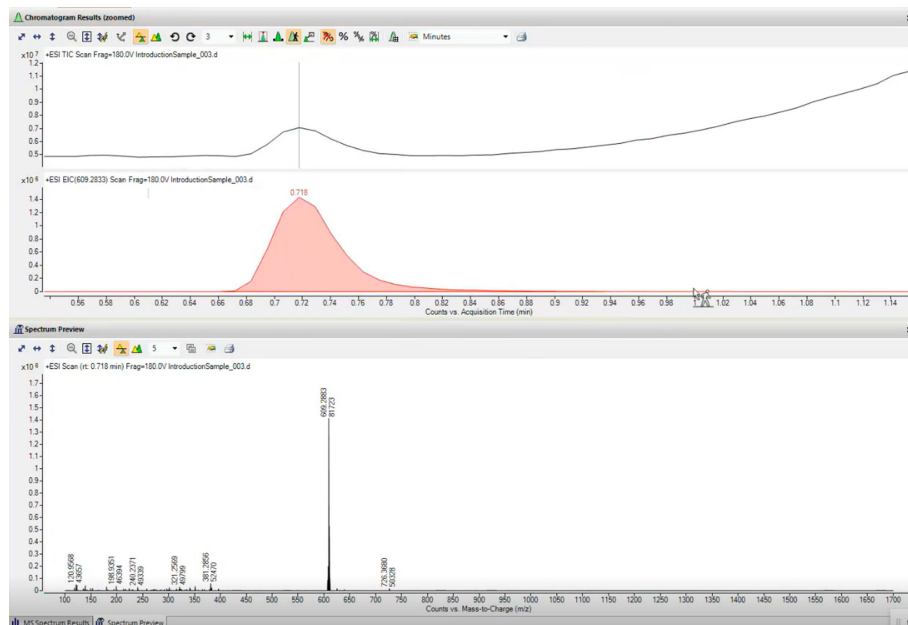
3. In the main window, in the Navigator View, right-click a chromatogram and select **Extract Chromatograms** from the shortcut menu.



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



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 Setting Up and Running a Sequence

A sequence can only be run after an acquisition method has been developed.

1. Click the **Sequence** layout.
2. In the Sequence window, enter the **Properties** information.
 - a. In the navigation pane at the left of the window, click **General > Properties**.
 - b. Select or clear **Audit Trail Activated**. Audit Trail cannot be turned off once it has been enabled.
 - c. Enter the **Description** for the sequence.

The screenshot shows the 'Sequence Properties' dialog box with the 'General > Properties' tab selected. The left navigation pane shows 'General' expanded, with 'Properties' selected. The main area contains a 'Location' field, a 'Version' field, a 'Show Audit Trail' button, and an 'Audit Trail Activated' checkbox. Below these is a large 'Description' text area. At the bottom, there are fields for 'Created at', 'Modified at', 'Created by' (SYSTEM (SYSTEM)), and 'Modified by'.

3. Enter the **Run Options** information:
 - a. In the navigation pane at the left of the window, click **General > Run Options**.
 - b. Select or clear **Use Barcode Reader Before Injection**.

- c. Click either **Inject anyway** or **Abort current injection**.

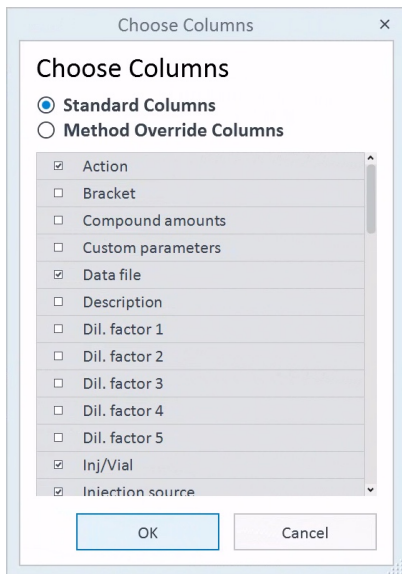
The screenshot shows the 'Run Options' dialog box. On the left is a navigation pane with three items: 'General Properties', 'Run Options' (which is highlighted in blue), and 'Injections Table'. The main area of the dialog is titled 'Injection report printer:' with a dropdown menu showing '[Printer specified in processing method]'. Below this is a section titled 'Barcode Reader' which contains a checkbox 'Use Barcode Reader Before Injection' and a sub-section 'On Barcode Mismatch' with two radio buttons: 'Inject anyway' (which is selected) and 'Abort current injection'. At the bottom is a section titled 'Data processing error handling for High Resolution Mass Spectra Data' which contains a checkbox 'Use Data processing error handling' and a sub-section 'On Data processing error' with two radio buttons: 'Ignore error and continue' (which is selected) and 'On error, pause after current injection'.

4. *Optional.* Add a column to the Sequence table.
- a. In the navigation pane at the left of the window, click **Injections > Table**.

Setting Up and Running a Sequence

Setting Up and Running a Sequence

- b. In the upper left corner of the Sequence table, click . The Choose Columns dialog opens.



- c. Select the desired column (or columns).
 - d. Click **Method Override Columns**.
 - e. Select the desired override column (or columns).
 - f. Click **OK**.
5. Enter the **Table** information:
 - a. In the navigation pane at the left of the window, click **Injections > Table**.
 - b. Enter the information for each injection.

	Action	Vial	Samp...	Run t...	Level	Acq...	Proc...	Inj/Vial	Volu...	Inject...	Samp...	Samp...	Data...	Label	Samp...
1	Inject	1	Sa...			default...		1	Use Me...	HipAls	0.00000				
2	Inject	1	Sa...			default...		1	Use Me...	HipAls	0.00000				

Setting Up and Running a Sequence

Setting Up and Running a Sequence

6. Select the rows to submit as part of the run. By default, all rows are selected.
7. To save your results to a sub-folder inside your Results folder, change the **Result Path** by clicking browse and creating a new folder. Results cannot be saved to an existing .rslt folder.

Result path	E:\CDSProjects\Learning Products\Results	>	...	Save result as	Separate single injections ▼
Result name		>			

8. The **Result name** defaults to the last filename used for this instrument and project. Enter a filename by typing in the field or using tokens. If left blank, the instrument name and local date & time is used for the filename.
9. Select to **Save result as**:
 - **One result set** to save the data acquired for all sample injections in the sequence within a single result set. For example:
 - ResultSetName: A1, B1, A2, B2, A3, B3, A4, B4, A5, B5
 - **Separate single injections** to save each sample injection in the sequence as an individual result set. When using this option, result set names will be based on the value of the **data file** column for each sequence row. For example:
 - DataFileName: A1
 - DataFileName: B1
 - DataFileName: A2
 - DataFileName: B2
10. *Optional.* Click Validate to validate the sequence against the specified acquisition method. The columns validated are **Acq. method**, **Volume**, and any override parameters. If invalid entries are found, correct those entries.
11. Click Run. If this is a new "Untitled" sequence, the sequence file is saved as the instrument name + the local date and time.
Submitted sequences are placed in the **Run Queue**, where runs can be stopped or paused, and sequences can be edited. A result folder (.rslt) is created when the run is added to the Run Queue.



7

Maintenance

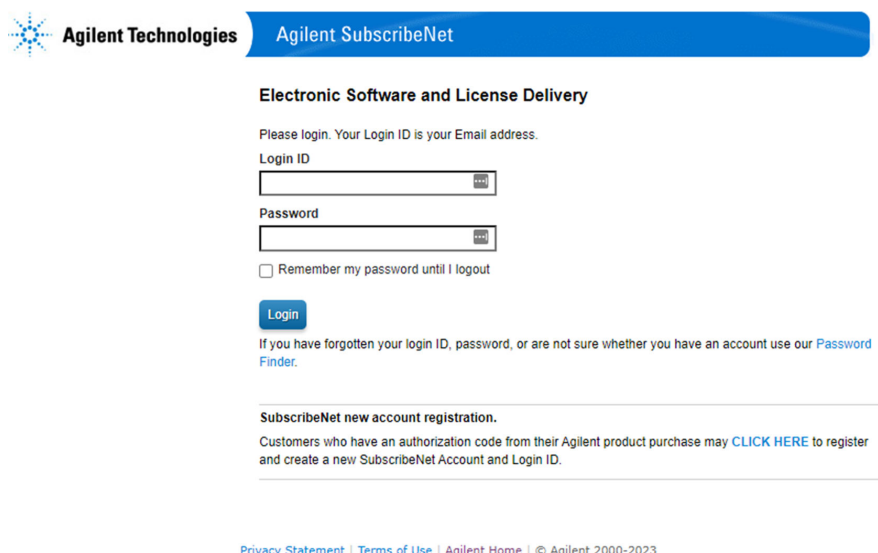
Register on Agilent SubscribeNet (New Account Registration)	62
Perform Back Up and Platform Best Practices	64

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.subscribe.net.com/>.



The screenshot shows the Agilent SubscribeNet website. At the top, there is a blue header with the Agilent logo and the text "Agilent Technologies" and "Agilent SubscribeNet". Below the header, the main content area is titled "Electronic Software and License Delivery". Under this title, there is a message: "Please login. Your Login ID is your Email address." Below this message are two input fields: "Login ID" and "Password". Below the "Password" field is a checkbox labeled "Remember my password until I logout". Below the checkbox is a blue "Login" button. Below the "Login" button is a link: "If you have forgotten your login ID, password, or are not sure whether you have an account use our [Password Finder](#)." Below this link is a section titled "SubscribeNet new account registration." followed by the text: "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." At the bottom of the page, there is a footer with links: "Privacy Statement | Terms of Use | Agilent Home" and copyright information: "© 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**Register on Agilent SubscribeNet (New Account Registration)**

- 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**.



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.

Maintenance

Perform Back Up and Platform Best Practices



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. 2026

G1960-90158

May 2026

