



Agilent 6475 Triple Quad LC/MS System

## User Guide



# Notices

## Document Identification

DocNo D0025685 Rev. A.01  
DE22440912  
April 2023

## Copyright

© Agilent Technologies, Inc. 2023

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

## Software Revision

This guide is valid for MassHunter 12.0, until superseded.

## Instrument Manufacturing

Manufactured by Agilent  
Technologies Singapore Pte. Ltd.  
No. 1 Yishun Avenue 7, Singapore  
768923



## Operating Temperature

Operating Temperature: 15°C to 35°C  
Storage Temperature: -40°C to 70°C

## 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.

# Content

<b>1</b>	<b>About this System</b>	<b>6</b>
	<b>Sustainability Information</b>	<b>7</b>
	Packaging End-of-Life	7
	Product End-of-Life	7
	<b>Agilent Regulatory Compliance Statement</b>	<b>8</b>
	CE Compliance	8
	UK Compliance	9
	Electromagnetic Compatibility	10
	EMC Declaration for South Korea	11
	Detachable Power Cord Declaration for Japan	11
	Sound Emission Certification for Federal Republic of Germany	11
	Waste Electrical and Electronic Equipment (WEEE) Directive	12
	<b>Technical Support</b>	<b>13</b>
	Agilent Community	13
<b>2</b>	<b>Hardware Overview</b>	<b>14</b>
	<b>6475A LC/TQ</b>	<b>15</b>
	Front view	15
	Calibrant Delivery System	16
	Side view	17
	<b>Instrument Status Indicator Lights</b>	<b>18</b>
	<b>Instrument Status Indicators in Acquisition</b>	<b>19</b>
	<b>Ion Sources</b>	<b>20</b>
	Agilent Jet Stream (AJS) source	21
	Electrospray Ionization (ESI) source	22
	Atmospheric Pressure Chemical Ionization (APCI) source	23
	Multimode Ionization (MMI) source	24
	Nanospray ESI source	25

## Content

<b>3</b>	<b>Tuning</b>	<b>26</b>
	<b>About Tuning</b>	<b>27</b>
	Autotune and Checktune	27
	<b>Starting the Instrument</b>	<b>28</b>
	<b>Calibrating the LC/TQ (Checktune)</b>	<b>29</b>
	<b>Calibrating the LC/TQ (Autotune)</b>	<b>30</b>
	<b>Generating a Detailed Tune Report</b>	<b>31</b>
	<b>Scheduling a Checktune</b>	<b>32</b>
	<b>Scheduling an Autotune</b>	<b>33</b>
<b>4</b>	<b>Basic Operation</b>	<b>34</b>
	<b>Starting Up and Shutting Down the Instrument</b>	<b>36</b>
	Starting in Standby mode	36
	Turning off the instrument	36
	<b>Starting the Acquisition Software</b>	<b>38</b>
	Acquisition Windows	39
	<b>Preparing the LC Modules</b>	<b>45</b>
	<b>Preparing the Triple Quad LC/MS</b>	<b>50</b>
	Performing a Checktune or Autotune	50
	Switching LC stream to MS	50
	Monitoring MS baseline and spectral displays	50
	Viewing the System Activity Log for events and errors	51

## Content

	<b>Setting Up an Acquisition Method</b>	<b>52</b>
	<b>Setting Up and Running a Single Sample</b>	<b>56</b>
	<b>Setting Up and Running a Worklist</b>	<b>57</b>
	<b>Reviewing Results with Qualitative Analysis</b>	<b>61</b>
	<b>Analyzing Data with Quantitative Analysis</b>	<b>62</b>
<b>5</b>	<b>Maintenance</b>	<b>63</b>
	<b>Routine Tasks</b>	<b>64</b>
	Calibrant Delivery System and Divert Valve	64
	Sources	65
	Nebulizer	68
	<b>Periodic Tasks</b>	<b>69</b>
	Calibrant Delivery System and Divert Valve	69
	Sources	70
	Nebulizer	75
	Corona Needle	80
	Ion Injector	82
	Rough Pump	85
	<b>Disinfection</b>	<b>90</b>



# 1

## About this System

Sustainability Information	7
Packaging End-of-Life	7
Product End-of-Life	7
Agilent Regulatory Compliance Statement	8
CE Compliance	8
UK Compliance	9
Electromagnetic Compatibility	10
EMC Declaration for South Korea	11
Detachable Power Cord Declaration for Japan	11
Sound Emission Certification for Federal Republic of Germany	11
Waste Electrical and Electronic Equipment (WEEE) Directive	12
Technical Support	13
Agilent Community	13

## Sustainability Information

### Packaging End-of-Life

The Agilent 6475 triple quadrupole LC/MS system is packaged using cardboard and a polyethylene bag. In the US, the cardboard is readily recyclable; while the plastic bag is recyclable, the infrastructure is not available in the majority of the market. In the US, Agilent is enrolled in the How2Recycle program to facilitate the communication of material recycling instructions. In select countries, Agilent participates in the EU/UK Green DOT Package Recycling Program to support the take-back and responsible management of the packaging materials at their end-of-life. In the remaining EU countries, the cardboard is readily recyclable while the plastic bag is recycled, landfilled, or incinerated depending on the country and its available infrastructure.

### Product End-of-Life

Agilent offers a take-back program for its 6475 triple quadrupole LC/MS system customers in US, EU, and UK markets.

# Agilent Regulatory Compliance Statement

## CE Compliance



Your Agilent instrument has been designed to comply with the requirements of the applicable directives of the European Union, such as Electromagnetic Compatibility (EMC) Directive, Low Voltage Directive (LVD), Machinery Directive (MD), RoHS Directive, etc.

Agilent has confirmed that each product complies with the relevant Directives by testing samples against the harmonized EN (European Norm) standards published on the Official Journal of the European Union (OJEU).

Proof that a product complies with these directives is indicated by:

- the CE Marking appearing on the rear of the product, and
- the documentation package that accompanies the product containing a copy of the Declaration of Conformity. The Declaration of Conformity is the legal declaration by Agilent that the product complies with the relevant directives listed above, and shows the EN standards to which the product was tested to demonstrate compliance.



## UK Compliance



Your Agilent instrument has been designed to comply with the requirements of the applicable regulations of the United Kingdom, such as The Electromagnetic Compatibility Regulations 2016, The Electrical Equipment (Safety) Regulations 2016, The Supply of Machinery (Safety) Regulations 2008, The Restriction of the Use of Certain Hazardous Substances in Electrical and Electronic Equipment Regulations 2012, etc.


Agilent has confirmed that each product complies with the relevant Regulations by testing samples against the designated standards published on GOV.UK.

Proof that a product complies with these regulations is indicated by:

- the UKCA Marking appearing on the rear of the product, and
- the documentation package that accompanies the product containing a copy of the Declaration of Conformity. The Declaration of Conformity is the legal declaration by Agilent that the product complies with the relevant regulations listed above, and shows the designated standards to which the product was tested to demonstrate compliance.

## Electromagnetic Compatibility

This product conforms to the following regulations on Electromagnetic Compatibility (EMC) and Radio Frequency Interference (RFI):

- CISPR 11/EN 55011: Group 1, Class A
- EC/EN 61326-1
- AUS/NZ 
- Canada ICES-001 (This ISM device complies with Canadian ICES-001. Cet appareil ISM est conforme à la norme NMB-001 du Canada).

**Group 1 ISM equipment** Group 1 contains all Industrial, Scientific and Medical (ISM) equipment in which there is intentionally generated and/or used conductively coupled radio- frequency energy which is necessary for the internal functioning of the equipment itself.

**Class A equipment** This equipment is suitable for use in all establishments other than domestic and those directly connected to a low voltage power supply network which supplies buildings used for domestic purposes.

This device complies with the requirements of CISPR11, Group 1, Class A as radiation professional equipment. Therefore, there may be potential difficulties in ensuring electromagnetic compatibility in other environments, due to conducted as well as radiated disturbances.

If this equipment does cause harmful interference to radio or television reception, which can be determined by turning the equipment off and on, the user is encouraged to try one or more of the following measures:

- 1 Relocate the radio or antenna.
- 2 Move the device away from the radio or television.
- 3 Plug the device into a different electrical outlet, so that the device and the radio or television are on separate electrical circuits.
- 4 Make sure that all peripheral devices are also certified.
- 5 Make sure that appropriate cables are used to connect the device to peripheral equipment.
- 6 Consult your equipment dealer, Agilent Technologies, or an experienced technician for assistance.

Changes or modifications not expressly approved by Agilent Technologies could void the user's authority to operate the equipment.

## EMC Declaration for South Korea

### 사용자안내문

This equipment has been evaluated for its suitability for use in a commercial environment. When used in a domestic environment, there is a risk of radio interference.

이 기기는 업무용 환경에서 사용할 목적으로 적합성평가를 받은 기기로서 가정용 환경에서 사용하는 경우 전파간섭의 우려가 있습니다 .

※ 사용자 안내문은 " 업무용 방송통신기자재 " 에만 적용한다 .

## Detachable Power Cord Declaration for Japan

電源コードセットの取扱いについて（日本国内向け）

製品には、同梱された電源コードセットをお使いください。同梱された電源コードセット

は、他の製品では使用できません。

Notice - The power cords for Japanese market

Your product must only use the power cord that was shipped with this product. Do not use this power cord with any other product.

## Sound Emission Certification for Federal Republic of Germany

### Sound pressure

Sound pressure  $L_p < 70$  dB(A) according to DIN EN ISO 7779.

### Schalldruckpegel

Schalldruckpegel  $LP < 70$  dB(A) nach DIN EN ISO 7779.

## Waste Electrical and Electronic Equipment (WEEE) Directive

This product complies with the European WEEE Directive marking requirements. The affixed label indicates that you must not discard this electrical/electronic product in domestic household waste.



### NOTE

Do not dispose of in domestic household waste.

To return unwanted products, contact your local Agilent office or see <https://www.agilent.com/environment/product/index.shtml> for more information.

## Technical Support

To find your local sales and support contact, visit the following page:

**<https://www.agilent.com/en/contact-us/page>**

For technical support, visit the following page:

**<https://www.agilent.com/en/support>**

## 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 new videos, documents, tools, and webinars relevant to your work.

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



## 2 Hardware Overview

6475A LC/TQ 15

Front view 15

Calibrant Delivery System 16

Side view 17

Instrument Status Indicator Lights 18

Instrument Status Indicators in Acquisition 19

Ion Sources 20

Agilent Jet Stream (AJS) source 21

Electrospray Ionization (ESI) source 22

Atmospheric Pressure Chemical Ionization (APCI) source 23

Multimode Ionization (MMI) source 24

Nanospray ESI source 25

# 6475A LC/TQ

## Front view

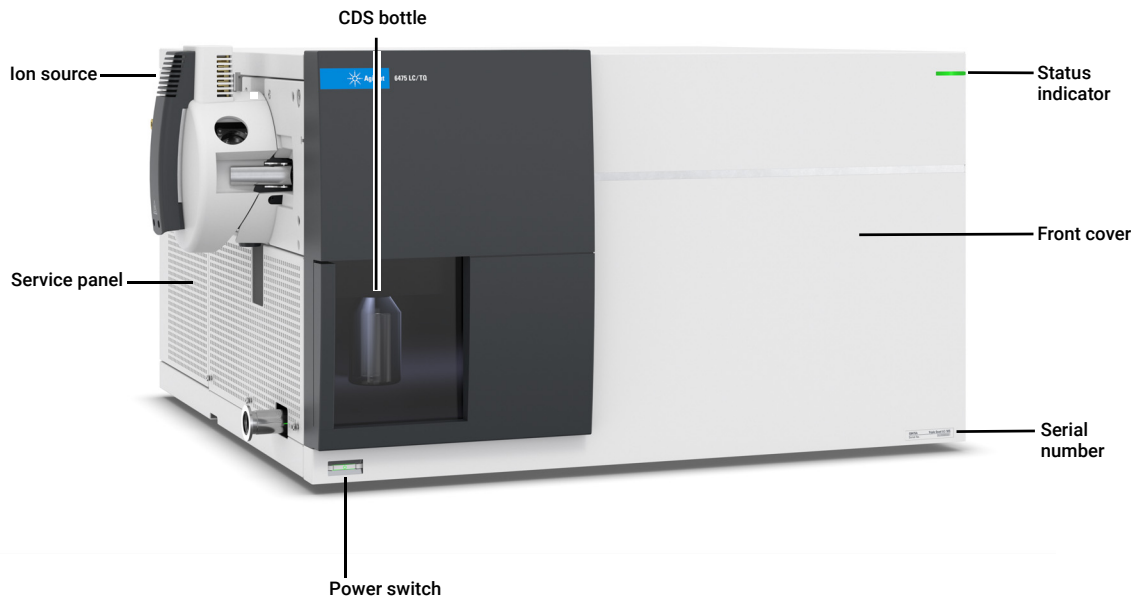


Figure 1. Front view of 6475A LC/TQ

## Calibrant Delivery System

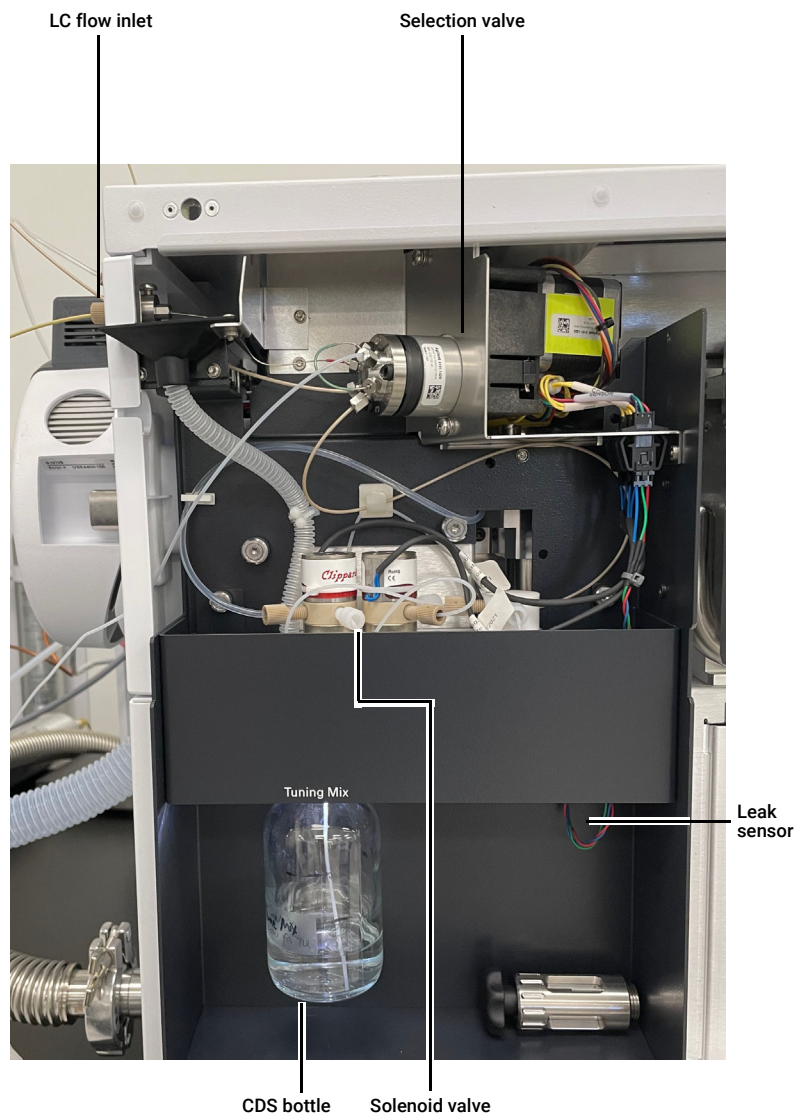


Figure 2. Calibrant delivery system



### Side view

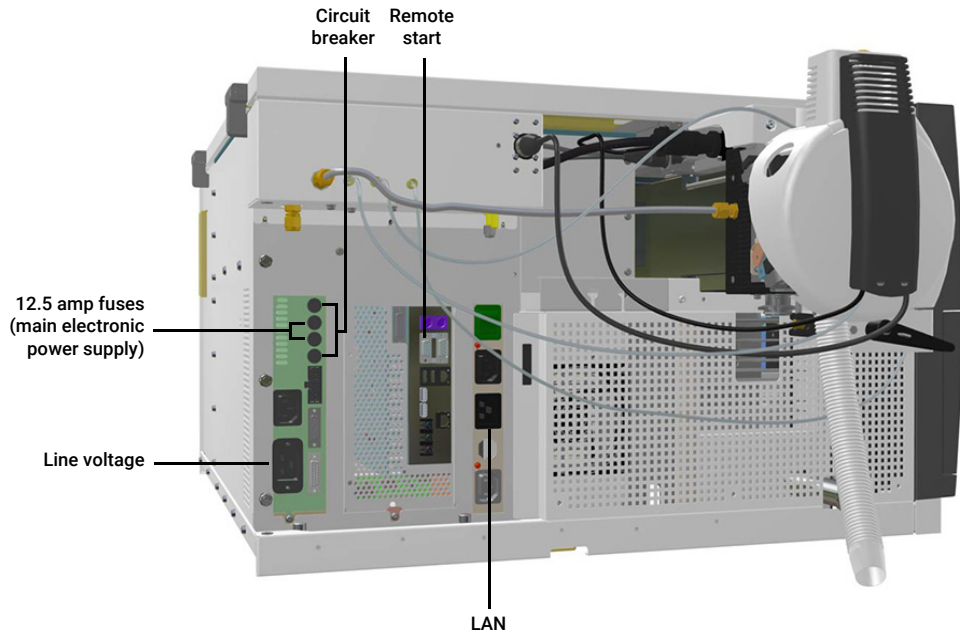


Figure 3. Side view of 6475A LC/TQ

# Instrument Status Indicator Lights

**Table 1** Status indicator lights on instrument

	<p>Instrument State = BOOTING (FPGA Loaded)            Status LED = Blinking Yellow            Power Switch LED = Yellow</p>		<p>Instrument State = ACQUISITION            Status LED = Blinking Green            Power Switch LED = Green</p>
	<p>Instrument State = BOOTING (LINUX Loaded)            Status LED = Yellow            Power Switch LED = Green</p>		<p>Instrument State = FAULT            Status LED = Red            Power Switch LED = Green</p>
	<p>Instrument State = PUMPING DOWN            Status LED = Blinking Yellow            Power Switch LED = Green</p>		<p>Instrument State = VENTING            Status LED = Blinking Red            Power Switch LED = Yellow</p>
	<p>Instrument State = STANDBY            Status LED = Yellow            Power Switch LED = Green</p>		<p>Instrument State = Vented            Status LED = Red            Power Switch LED = Yellow</p>
	<p>Instrument State = IDLE            Status LED = Yellow            Power Switch LED = Green</p>		<p>Instrument State = POWERING DOWN            Status LED = Off            Power Switch LED = Blinking Yellow</p>
	<p>Instrument State = BACKGROUND SCAN            Status LED = Green            Power Switch LED = Green</p>		<p>Instrument State = Off            Status LED = Off            Power Switch LED = Off</p>

## Instrument Status Indicators in Acquisition

In the Acquisition program, the Instrument Status window indicates the state of the instrument.

■ Gray - Offline: The instrument is configured with the system and available to use, but the AIC/workstation is not currently running. The amount of time for an instrument to reach the Idle state depends on each instrument.

■ Yellow - Unknown: The device is in an unknown run state.

■ Dark Yellow - Not Ready: The instrument is connected but is not ready to run (due to not reaching the correct temperature or pressure required by the method, for example).

■ Green - Idle: The instrument is on and ready to process samples.

■ Teal - Standby: The instrument is in a standby/sleep state. Put the instrument in Standby mode when it is not in use or when you want to change the ion source. When the instrument is in Standby mode:

- The source interface is on, the drying gas remains heated, and nebulizer flows are maintained.
- The source and ion optics voltages are turned off, and the mass spectrometer is not sending spectra to the computer.
- For APCI and multimode sources: Vaporizer gas heaters are turned down.
- For the Agilent Jet Stream (AJS) source: The sheath gas temperature, drying gas flow, and drying gas temperature are reduced.
- The spray chamber high voltages are turned off.
- The mass spectrometer stops generating spectra.

■ Purple - Pre-run/Injecting: The instrument is on and is preparing to start acquisition.

■ Blue - Running, Post Run: The instrument is currently collecting data.

■ Orange - Tuning: The instrument is in the process of adjusting MS parameters.

■ Light Orange - Not Ready: The device is not ready.

■ Red - Instrument Error: The instrument has an error and cannot process samples.

Blank - Disconnected: The connection to the instrument has been closed.

## Ion Sources

The following are supported ion sources:

- **“Agilent Jet Stream (AJS) source”** on page 21
- **“Electrospray Ionization (ESI) source”** on page 22
- **“Atmospheric Pressure Chemical Ionization (APCI) source”** on page 23
- **“Multimode Ionization (MMI) source”** on page 24
- **“Nanospray ESI source”** on page 25

For ion source maintenance, see **“Sources”** on page 65.

## Agilent Jet Stream (AJS) source

The Agilent Jet Stream (AJS) source uses the same ionization technique as the Electrospray source.

The AJS source uses Thermal Gradient Focusing Technology, which is a process in which super-heated nitrogen ( $N_2$ ) is used to improve ion generation and desolvation.

AJS source improves sensitivity 5x or more for many small-molecule compounds that undergo electrospray ionization.



**Figure 4.** AJS cover, ion source, and spray shield

## Electrospray Ionization (ESI) source

Electrospray relies in part on chemistry to generate analyte ions in solution before the analyte reaches the mass spectrometer. The LC eluent is sprayed (nebulized) into a chamber at atmospheric pressure in the presence of a strong electrostatic field and heated drying gas.

The electrostatic field causes further dissociation of the analyte molecules. The heated drying gas causes the solvent in the droplets to evaporate. As the droplets shrink, the charge concentration in the droplets increases. Eventually, the repulsive force between ions with like charges exceeds the cohesive forces and ions are ejected (desorbed) into the gas phase. These ions are attracted to and pass through a capillary sampling orifice into the mass analyzer.



**Figure 5.** ESI cover, ion source, and spray shield

## Atmospheric Pressure Chemical Ionization (APCI) source

In Atmospheric Pressure Chemical Ionization (APCI), the LC mobile phase is sprayed through a heated vaporizer (typically 250°C to 400°C) at atmospheric pressure. The heat vaporizes the liquid. The resulting gas-phase solvent molecules are ionized by electrons discharged from a corona needle.

The solvent ions then transfer charge to the analyte molecules through chemical reactions (chemical ionization). The analyte ions pass through a capillary sampling orifice into the mass analyzer. APCI is applicable to a wide range of polar and nonpolar molecules.

APCI rarely results in multiple charging, so APCI is typically used for molecules smaller than 1,500 u.

Because high temperatures are also involved, APCI is not appropriate for analysis of large biomolecules. APCI is used with normal-phase chromatography more often than electrospray is because the analytes are usually nonpolar.

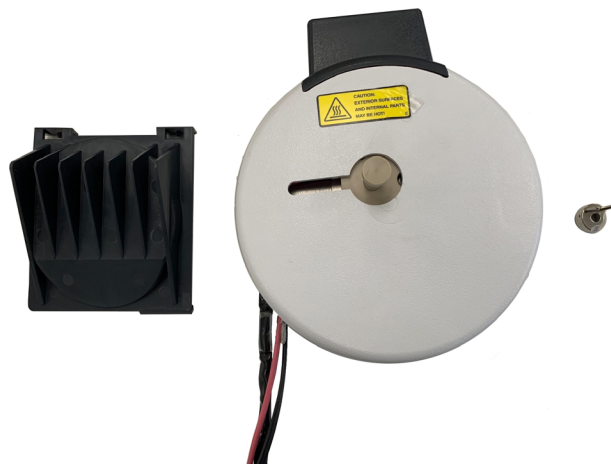


**Figure 6.** APCI cover, ion source, and spray shield

## Multimode Ionization (MMI) source

The Multimode source for LC/MS can simultaneously do ESI and APCI ionization. The Multimode source can operate in ESI mode only, in APCI mode only, or in mixed ESI/APCI mode. This technology significantly improves the speed, accuracy, and productivity of high-throughput screening in drug discovery and other research applications.

High-throughput screening is used to analyze large numbers of chemical compounds rapidly and with high sensitivity. In order to be analyzed by mass spectrometry, all compounds must be converted into ions. Since not all molecules respond to the same mode of ionization, multiple ionization sources were developed to facilitate high-throughput screening.



**Figure 7.** MMI cover, ion source, and spray shield

Simultaneous ESI and APCI capability and high LC flow rate compatibility (up to 2 mL/min) means more compounds can be detected in less time with the Agilent Multimode source.



## Nanospray ESI source

The Agilent nanospray ESI source is a flexible solution for laboratories running low-flow LC/MS applications. The nanospray source allows both commercially available and custom nanoflow columns to be used with Agilent LC/MS systems.

The spray position can be configured in orthogonal, direct, or direct off-axis orientations, to best suit your application. Visualize your spray stability with built-in video capture displayed on the control PC monitor. Reference mass solution can also be introduced to ensure the most accurate performance on the Agilent TOF and Q-TOF instruments.



**Figure 8.** Nanospray ESI source



## 3

# Tuning

About Tuning 27

Autotune and Checktune 27

Starting the Instrument 28

Calibrating the LC/TQ (Checktune) 29

Calibrating the LC/TQ (Autotune) 30

Generating a Detailed Tune Report 31

Scheduling a Checktune 32

Scheduling an Autotune 33

## About Tuning

### Autotune and Checktune

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 verifies peak width and mass axis to make sure they are correct before you start your acquisition. This checktune takes approximately 3 minutes to run in each polarity. See **“Calibrating the LC/TQ (Checktune)”** on page 29.

**Autotune** only needs to run after preventative maintenance or if you find a problem with Checktune. An Autotune can take approximately 15 to 20 minutes for each polarity. Everything is automatic since the tuning mix is delivered by the calibrant delivery system, which is switched on automatically during the tune. See **“Calibrating the LC/TQ (Autotune)”** on page 30.

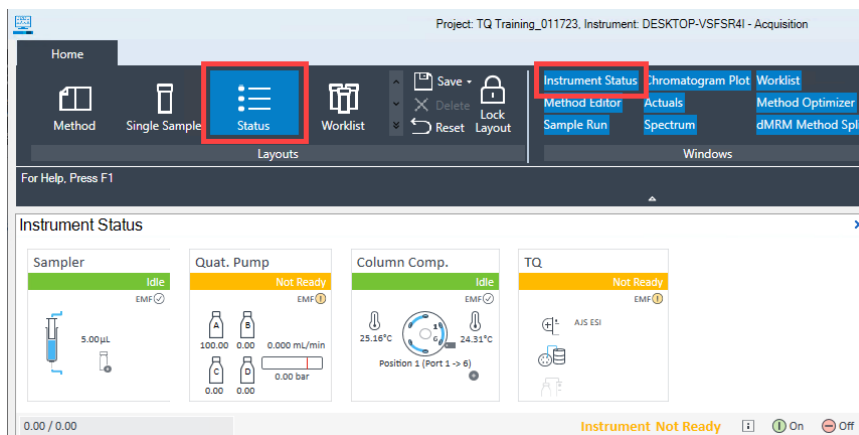
Use **Table 2** to select the correct tuning mix for the source installed.

**Table 2** Tuning mixes

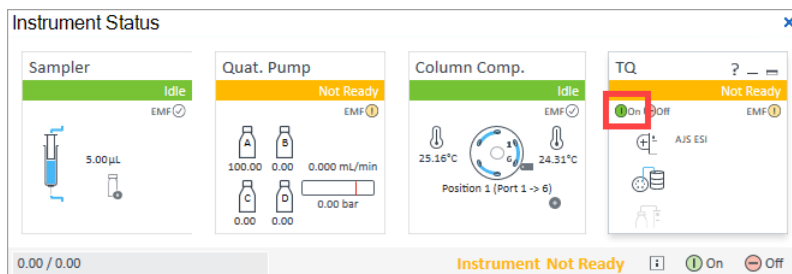
Source	Tuning mix
ESI	ESI-L Low Concentration Tuning Mix
AJS ESI	ESI-L Low Concentration Tuning Mix
MMI	MMI-L Low Concentration Tuning Mix
APCI	APCI-L Low Concentration Tuning Mix

## Starting the Instrument

- 1 Follow the procedures in **“Starting Up and Shutting Down the Instrument”** on page 36 to begin operation of the instrument.
- 2 Display the Instrument Status window if it is not visible in the Acquisition program, in one of the following ways:
  - Open the Status layout.
  - On the ribbon, click **Instrument Status** in the **Windows** section.




- 3 Hover over the TQ device pane and click **On**.

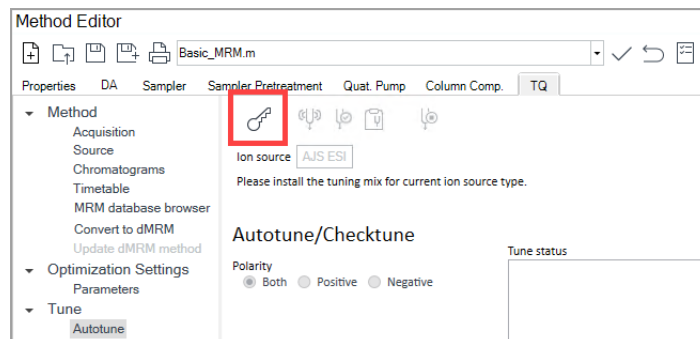


- 4 Monitor baseline and adjust plot if necessary. See **“Monitoring baseline and adjusting plot”** on page 41.

## Calibrating the LC/TQ (Checktune)



A Checktune can be run with the following ion sources: ESI, AJS ESI, MMI, and APCI.

- 1 In the Method Editor window, click the **TQ** tab.
- 2 Click **Tune > Autotune** in the left pane.
- 3 Click  **Request tune control**.




This button locks control of the TQ instrument. You cannot start a single sample run or a worklist when Tune has control of the TQ instrument.

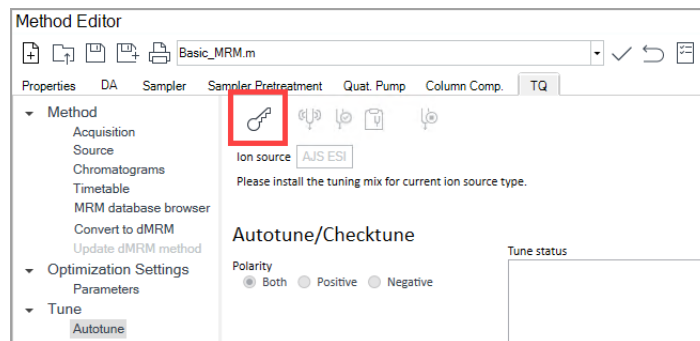
When clicked, tune setpoints are applied. If you do not select an action within the timeout period, the system will go back to the method setpoints.

- 4 Select a polarity. If **Both** is selected, the system will automatically run in the positive mode followed by the negative mode.
- 5 Click  **Checktune the instrument**.
- 6 When the tune completes, review the report.
- 7 Click  **Release tune control** to release control of the TQ instrument.

## Calibrating the LC/TQ (Autotune)



An Autotune can be run with the following ion sources: ESI, AJS ESI, MMI, and APCI.

- 1 In the Method Editor window, click the **TQ** tab.
- 2 Click **Tune > Autotune** in the left pane.
- 3 Click  **Request tune control**.




This button locks control of the LC/TQ instrument. You cannot start a single sample run or a worklist when Tune has control of the LC/TQ instrument.

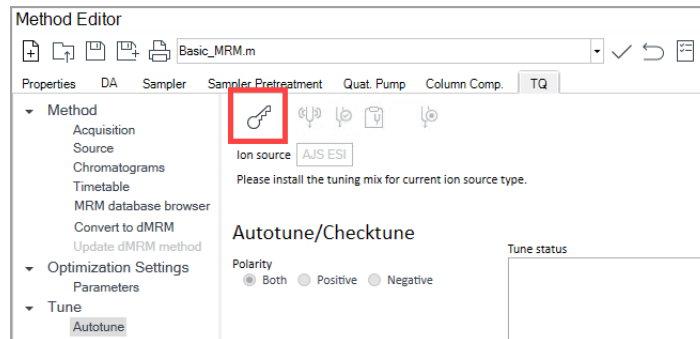
When clicked, tune setpoints are applied. If you do not select an action within the timeout period, the system will go back to the method setpoints.

- 4 Select a polarity.
  - If only **Positive** or **Negative** polarity is selected, subsequence acquisition methods are only available with that polarity.
  - If **Both** is selected, the system will automatically run in the positive mode followed by the negative mode.
- 5 Click  **Autotune the instrument**.
- 6 When the tune completes, review the report.
- 7 Click  **Release tune control** to release control of the LC/TQ instrument.



# Generating a Detailed Tune Report

A detailed tune report can be generated after you have run Autotune or Checktune.

- 1 In the Method Editor window, click the **TQ** tab.
- 2 Click **Tune > Autotune** in the left pane.
- 3 Click  **Request tune control**.



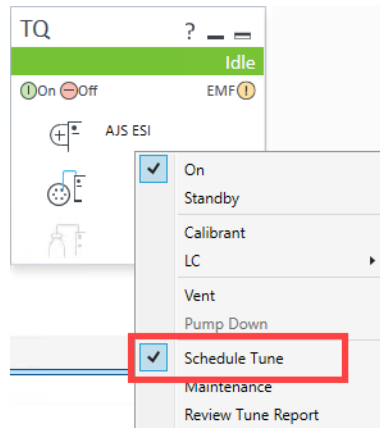
This button locks control of the TQ instrument. You cannot start a single sample run or a worklist when Tune has control of the TQ instrument.

- 4 Click  **Generate Detailed Tune Report**. If this button is not available, then you need to run either a Checktune or an Autotune first.
- 5 Click  **Release tune control** in the toolbar to release control of the TQ instrument.

## Scheduling a Checktune

To specify to run a Checktune automatically,

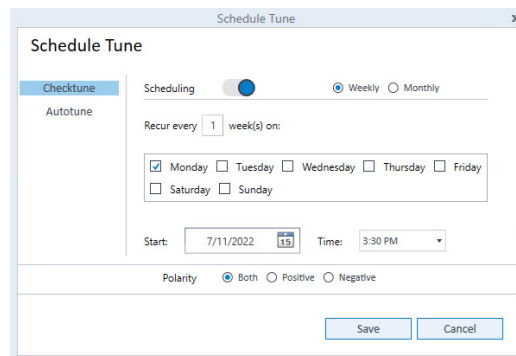
- 1 In the Instrument Status window, right-click the TQ device.
- 2 Click **Schedule Tune**.



The **Schedule Tune** dialog box opens.

- 3 Select **Checktune** in the left pane.

The right pane shows the information for scheduling a Checktune.



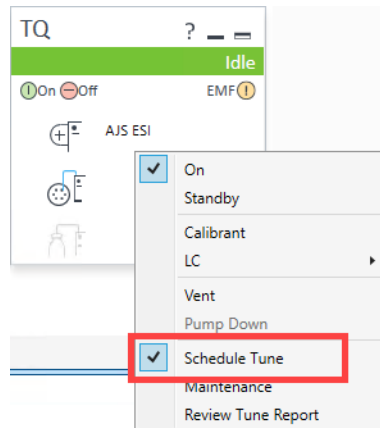
- 4 Click the **Scheduling** slider to switch **Scheduling** on.
- 5 Select either **Weekly** or **Monthly**.
- 6 Select other options to indicate how often to schedule the Checktune.
- 7 Click **Save**.



## Scheduling an Autotune

To specify to run an Autotune automatically,

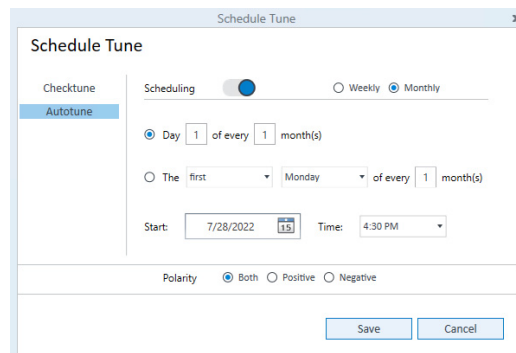
- 1 In the Instrument Status window, right-click the TQ device.
- 2 Click **Schedule Tune**.



The Schedule Tune Dialog Box opens.

- 3 Select **Autotune** in the left pane.

The right pane shows the information for scheduling an Autotune.



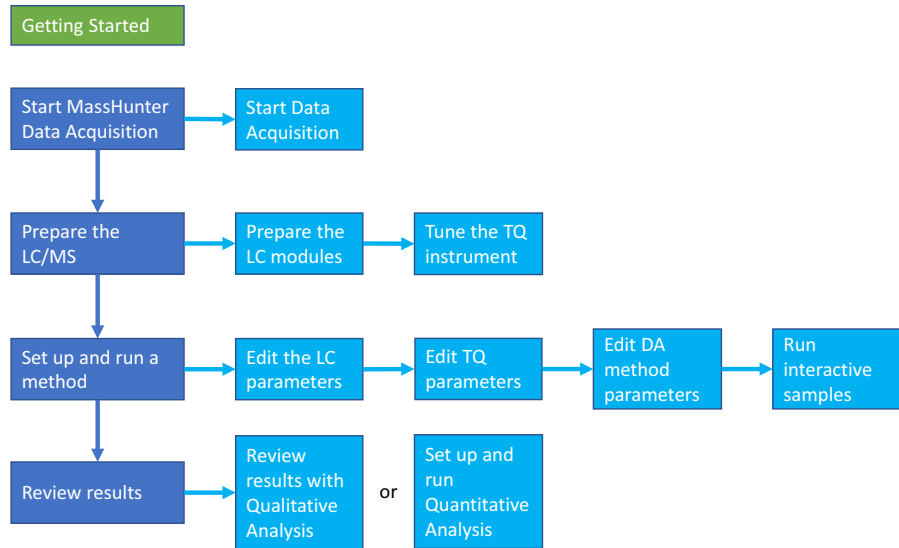
- 4 Click the **Scheduling** slider to switch **Scheduling** on.
- 5 Select either **Weekly** or **Monthly**.
- 6 Select other options to indicate how often to schedule the Autotune.
- 7 Click **Save**.

## 4

# Basic Operation

Starting Up and Shutting Down the Instrument	36
Starting in Standby mode	36
Turning off the instrument	36
Starting the Acquisition Software	38
Acquisition Windows	39
Preparing the LC Modules	45
Preparing the Triple Quad LC/MS	50
Performing a Checktune or Autotune	50
Switching LC stream to MS	50
Monitoring MS baseline and spectral displays	50
Viewing the System Activity Log for events and errors	51
Setting Up an Acquisition Method	52
Setting Up and Running a Single Sample	56
Setting Up and Running a Worklist	57
Reviewing Results with Qualitative Analysis	61
Analyzing Data with Quantitative Analysis	62

The roadmap below shows you the steps to set up and run a batch of samples from start to finish. Follow the instructions in this section to get started and to learn where to find the information to help you with each step in this roadmap.



## Starting Up and Shutting Down the Instrument

### Starting in Standby mode

- 1 Check that nitrogen gas for the drying gas and for the collision cell are turned on.
- 2 Turn on the power for the computer and monitor.
- 3 Check that the instrument and rough pump are plugged into a power outlet.
- 4 Press the front power switch on the instrument.

The vacuum system automatically starts to pump down the instrument, and the electronics are turned on. The system card boots up and the firmware starts.

### Turning off the instrument

Before you turn off the instrument, complete any acquisition steps and save your data.

#### CAUTION

**To prevent instrument damage, always vent the system before pressing the power button.**

#### WARNING

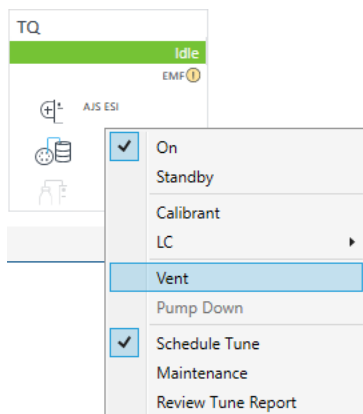
**Do not touch the spray shield and related spray chamber components. They are likely to be hot.**

- 1 Turn off the solvent flow.  
If you have any analyte in any of the components in the sample delivery system, flush the delivery system with pure solvent before you continue.
- 2 Check that the ballast valve on the rough pump is closed (if applicable).

## Basic Operation

### Turning off the instrument

- In Acquisition, right-click the TQ icon in the Instrument Status window and select **Vent** to vent the system. Click **Yes** to continue.



When you vent the system, the following occurs:

- The spray chamber high voltages, the drying gas heater, the nebulization flow, the detector, and other lens voltages are turned off.
  - The drying gas flow is set to Standby.
  - If the APCI source is installed, the vaporizer heater is also turned off.
  - If the installed source is an Agilent Jet Stream source, the sheath gas heater is also turned off.
  - The Rough Pump turns off when the turbo speed is below 20 percent.
- Wait until the instrument completely vents (reaching 760 Torr).
  - Close the Acquisition program and shut down the computer.
  - Press the power switch located in the lower left front corner of the instrument.



- Unplug the main power cable behind the service panel. This action prevents the turbo pump components from overheating.

## Starting the Acquisition Software

The instructions below include the following assumptions:

- The hardware and software are installed.
- The instrument is configured. For information on configuring your instrument, refer to the MassHunter Control Panel online Help.
- The LC modules and the Triple Quad LC/MS are turned on, but the LC pump is not running.

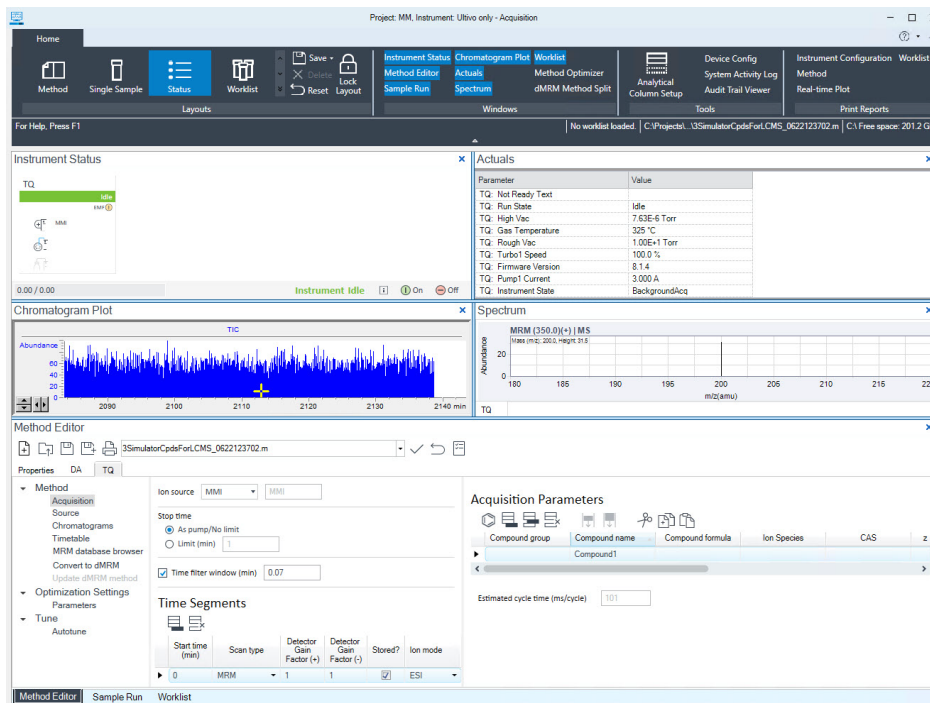
After installation, all the Agilent MassHunter Workstation Software icons appear on your desktop.

To start the Acquisition program,

- 1** Double-click the **Control Panel** icon on your desktop, or select **Windows Start > Agilent Technologies > Control Panel**.
- 2** In Control Panel, select an instrument and click **Launch** to open the project in Acquisition. For more information, refer to the MassHunter Control Panel online Help.

## Acquisition Windows

When the Acquisition program opens, the main window appears. Almost all of your work is done within the windows of this main window. These windows provide the tools to set up acquisition methods, run samples interactively or automatically, monitor instrument status, monitor runs, and tune the instrument.



### Showing and hiding the windows

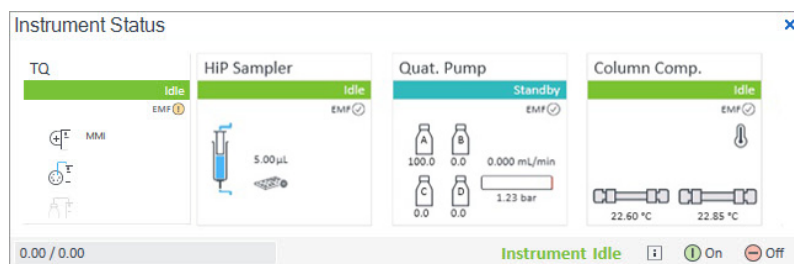
To show or hide a window, click the window name in the Windows group of the ribbon. Or, hide a window by clicking the **X** icon in the upper right corner of the window. At least one window must be shown.

Click in a window and Press **F1** to obtain help on the active window. Drag a window border to resize the window. If you double-click the title of the window, the window "floats" outside of the main window. Double-click the title bar again to "dock" the window.

#### Instrument Status window

This window displays the status of each device configured with the instrument: **Error**, **Not ready**, **Pre-run**, **Post-run**, **Running**, **Injecting**, **Idle**, **Offline**, or **Standby**. You also set non-method control and configuration parameters for the LC devices and the MS instrument.

This window displays the current status of each device both as text and by its color-coding. See **“Instrument Status Indicators in Acquisition”** on page 19.



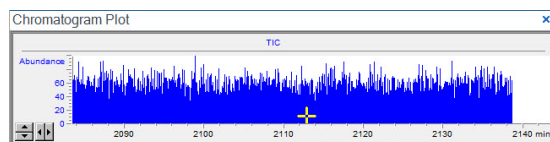
#### Actuals window

This window displays the current value of selected instrument parameters. See **“Set up to view real-time parameter values (actuals).”** on page 46 for more information.

Parameter	Value
TQ: Not Ready Text	
TQ: Run State	Idle
TQ: High Vac	7.63E-6 Torr
TQ: Gas Temperature	325 °C
TQ: Rough Vac	1.00E+1 Torr
TQ: Turbo1 Speed	100.0 %
TQ: Firmware Version	8.1.4
TQ: Pump1 Current	3.000 A
TQ: Instrument State	BackgroundAcq

#### Chromatogram Plot window

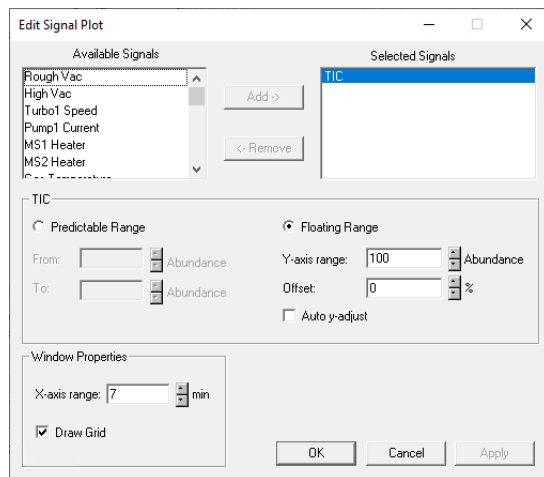
This window displays the chromatogram plots in real time. These plots can be user-defined signals and/or instrument parameters. You select the plots in the Chromatogram tab in the **QQQ** tab in the Method Editor window.





**Monitoring baseline and adjusting plot** The **Edit Signal Plot** dialog allows you to select various instrument parameters and signals for display in the Chromatogram Plot window. The signals and their display characteristics can be adjusted as described in this section. The customizations you make to the plot remain, even if you close the Acquisition software and reopen it.

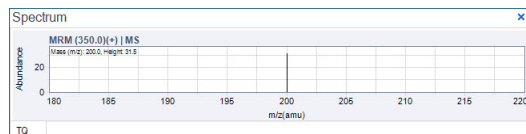
- 1 In the Chromatogram Plot window, right-click the Chromatogram Plot window, and select **Change**. The **Edit Signal Plot** dialog box opens.



- 2 Select the signal to monitor and set the X and Y range for the plot.
- 3 To add other MS signals, such as EICs, setpoints, and actuals, select these on the **TQ** tab in the Method Editor window.
- 4 Click **OK** to close the **Edit Signal Plot** dialog box.

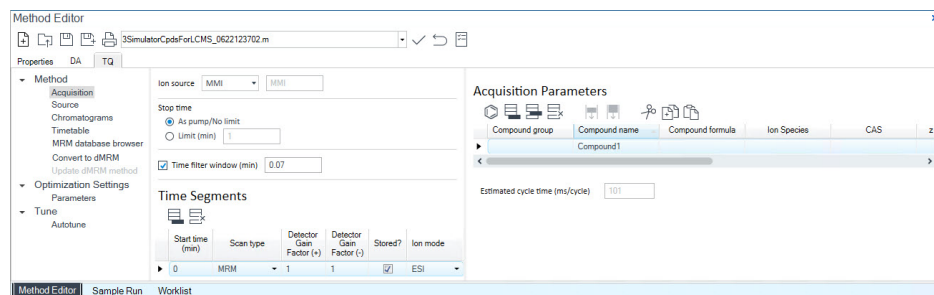
## Spectrum window

This window displays the spectral plot in real time. A different tab is available for different devices.



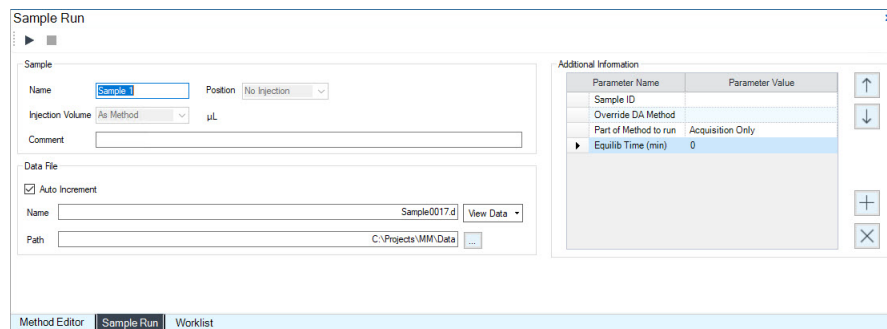
#### Method Editor window

In this window, enter acquisition parameters for the method. Tune values are displayed on the Autotune tab.



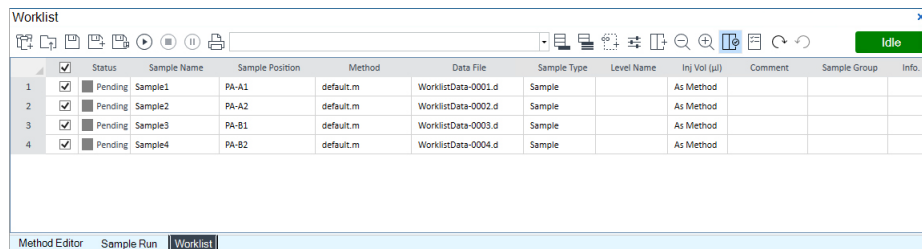
#### Sample Run window

In this window, enter sample information to run individual samples interactively, and start a single sample run. Specify an **Override DA Method** and select either **Both Acquisition and DA** (to run Data Analysis as part of the method) or **Acquisition Only** for the **Part of Method to run**.



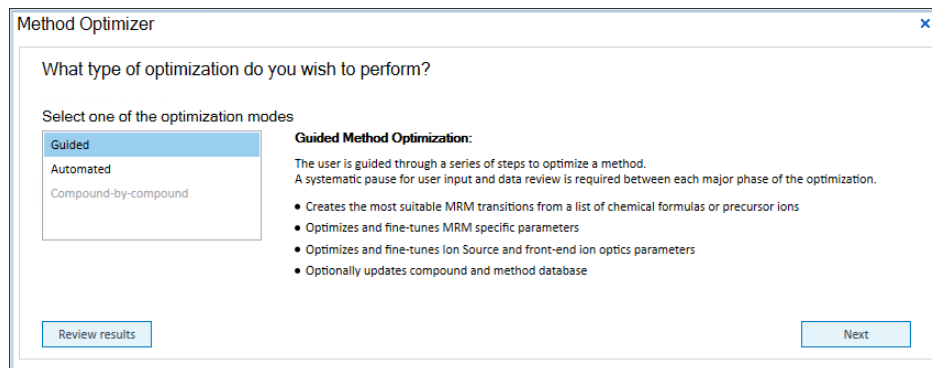
### Worklist window

In this window, enter sample information for multiple samples. When you run the worklist, the samples are automatically run in the order listed in the worklist. Select whether to run **Acquisition Only** or **Both Acquisition and DA** for the **Part of method to run** option in the Worklist Run Parameters dialog box.



### Method Optimizer window

Method Optimizer operates in both Workstation and Networked Workstation configurations. The Method Optimizer can automate MRM product ion selection, collision energies, fragmentor voltages, and source parameters.



**dMRM Method Split window**

The dMRM Method Split window operates in both Workstation and Networked Workstation configurations. This window facilitates splitting methods with many dMRM transitions into separate methods with fewer dMRM transitions.

dMRM Method Split

Please choose a method to split

...

Number of methods  Split

Maximum concurrent MRMs

Minimum dwell time (ms)

## Preparing the LC Modules

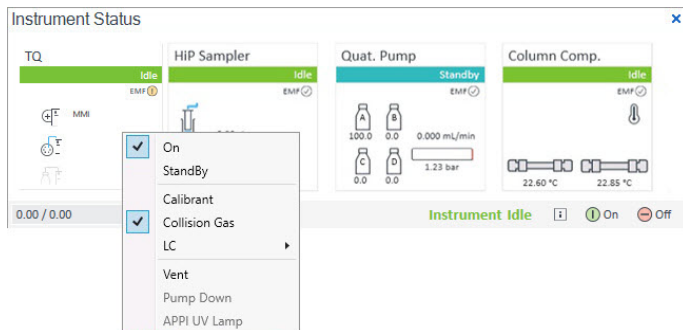
Read and follow the instructions in the online Help for each of the tasks in the checklist described on the following pages. While you condition or equilibrate the column, you can tune the Triple Quad LC/MS.

### 1 Switch LC stream to **Waste**.

When you are not acquiring data, you switch the direction of the LC stream away from the MS ion source and to waste.

If you have the LC connected to a VWD or DAD, you can still monitor the fluctuations of the VWD or DAD real-time chromatogram before a run.

- a Right-click the **TQ** device in the Instrument Status window.
- b Select **LC > Waste**.

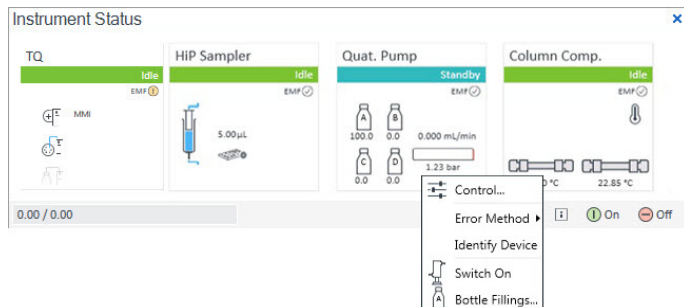


- 2 Purge the LC pump. Follow the directions for purging the pump in the *User Guide* for your pump.
- 3 Set up to condition or equilibrate the column.
  - a In the Method Editor window, click **Apply** in the toolbar to download the LC parameters to the LC.

## Basic Operation

### Preparing the LC Modules

- b Right-click an LC module in the Instrument Status window and select one of the commands to change any non-method control parameters, if needed.



- c Monitor the baseline and adjust the plot to make sure the column is equilibrated and the baseline stable. (See **step 4** and **step 5** in this procedure.)
- 4 Set up to view real-time parameter values (actuals).

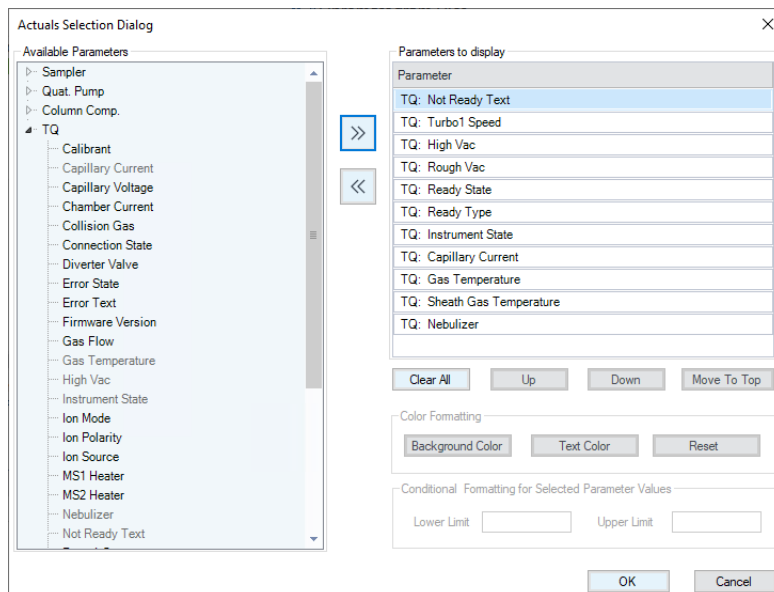
The Actuals window can be customized to display instrument conditions that you may want to view.

- a Click **Actuals** in the ribbon to display the Actuals window.
- b Right-click the **Actuals** window and select **Setup**.

Parameter	Value
TQ: Not Ready Text	
TQ: Run State	
TQ: High Vac	7.63E-6 Torr
TQ: Gas Temperature	325 °C
TQ: Rough Vac	1.00E+1 Torr
TQ: Turbo1 Speed	100.0 %
TQ: Firmware Version	8.1.4
TQ: Pump1 Current	3.000 A
TQ: Instrument State	BackgroundAcq

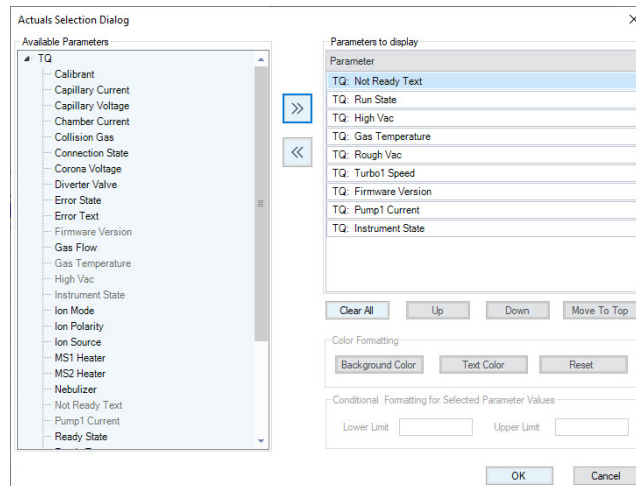
- c Select parameters to appear in the Actuals window.

In the **Available Parameters** list, expand a device to see a list of the instrument parameters for that device. Only configured devices will be shown in the list.



- d Select the parameters of interest, and click **>> Add**. Repeat this step for all the parameters you want to view.
- e Select any parameter in the **Parameters to display** list that you do not want included, and click **<< Remove**.
- f Customize the background color and the text color for any parameters in the **Parameters to display** list.
- g Set a conditional formatting range for the parameters that are numbers.

If the value of the parameter is not within the limits you entered, then the background of the parameter is set to red.



- h** Click **OK** when you finish selecting parameters.

The parameters you selected appear in the Actuals window. Note that the customizations you have made will remain, even if you close the Acquisition software and reopen it.



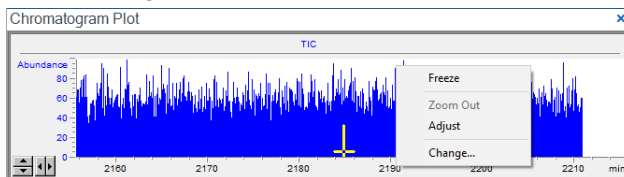
## Basic Operation

### Preparing the LC Modules

#### 5 Set up real-time plot displays.

As you condition the column, you can set up the displays to monitor the effluent.

- a In the Chromatogram Plots window, right-click the chromatogram plot, and select **Change**.



- b In the **Edit Signal Plot** dialog box, select the type of display signal, and click **OK**.

The screenshot shows the "Edit Signal Plot" dialog box. It has two panes: "Available Signals" and "Selected Signals". The "Available Signals" list includes Rough Vac, High Vac, Turbo1 Speed, Pump1 Current, MS1 Heater, and MS2 Heater. The "Selected Signals" list contains TIC. Below the panes, there are controls for the "TIC" signal. The "Predictable Range" radio button is unselected, and the "Floating Range" radio button is selected. The "From:" and "To:" fields are empty, and the "Y-axis range:" is set to 100 Abundance. The "Offset:" is set to 0%. There is an "Auto y-adjust" checkbox which is unchecked. At the bottom, there are "DK", "Cancel", and "Apply" buttons.

## Preparing the Triple Quad LC/MS

### Performing a Checktune or Autotune

See **Chapter 3**, “Tuning”.

### Switching LC stream to MS

After you condition the column and tune the instrument, switch the LC stream from **Waste** to **MS**. See “**Switch LC stream to Waste.**” on page 45 for how to do this.

### Monitoring MS baseline and spectral displays

If you did not monitor the LC baseline with a VWD or DAD, make sure that the instrument baseline is stable and no spectra of interfering intensity appear in the display.

If you did monitor the LC baseline with a VWD or DAD, change back to the default instrument displays.

- 1 Right-click the chromatogram plot, and click **Change**.
- 2 Select the MS signal, and click **OK**.

## Viewing the System Activity Log for events and errors

The System Activity Log provides information regarding errors to assist with troubleshooting.


To view the log, click **System Activity Log** in the ribbon of the Acquisition program, and view the logged events.

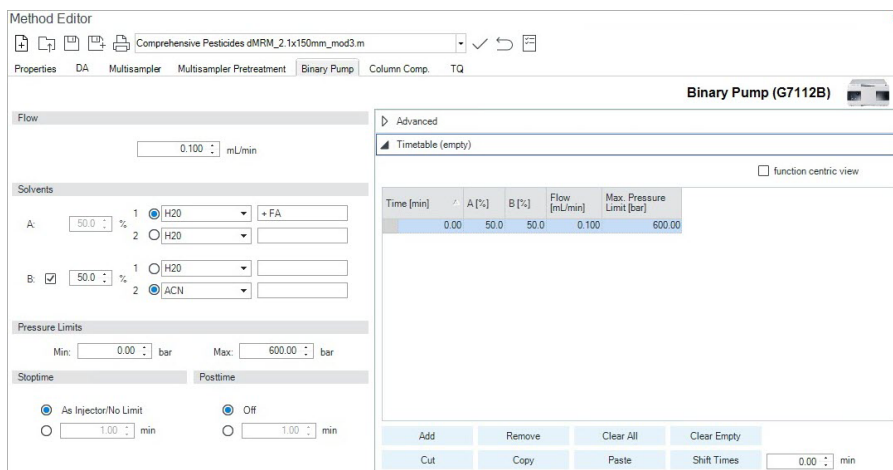


Date/Time	User	Description
2022-12-14 12:27:49-08:00	SYSTEM (SYSTEM)	Activity Log Viewer: User 'SYSTEM' has logged in.
2022-12-14 12:24:53-08:00	SYSTEM (SYSTEM)	Acquisition Engine: Method 'D:\Projects\GMTest\Methods\CompoundOptm (Version: NA)' loaded on instrument.
2022-12-14 12:24:47-08:00	SYSTEM (SYSTEM)	Acq Console: User has loaded layout 'Method (sys)'.
2022-12-14 12:24:33-08:00	SYSTEM (SYSTEM)	Acq Console: User 'SYSTEM' has logged in.
2022-12-14 12:24:31-08:00	SYSTEM (SYSTEM)	Engine Launcher: System Engines started.
2022-12-14 12:23:56-08:00	SYSTEM (SYSTEM)	Activity Log Enabled
2022-12-14 12:23:51-08:00	SYSTEM (SYSTEM)	User "SYSTEM (SYSTEM)" logged in
2022-12-13 17:39:36-08:00		OLSS host windows services started
2022-12-13 17:37:32-08:00		OLSS host windows services stopped
2022-12-08 11:27:20-08:00	SYSTEM (SYSTEM)	Activity Log Enabled
2022-12-08 11:27:17-08:00	SYSTEM (SYSTEM)	User "SYSTEM (SYSTEM)" logged in

1 - 11 of 11

## Setting Up an Acquisition Method

- 1 Set up the method in the Method Editor window:
  - a Enter or select the values and settings for each of the tabs below.
  - b *Optional.* To download the settings to the instrument, click ✓ **Apply**.
  - c To save the method, click  **Save As**.
  - d Name the method and click **OK**.
- 2 Enter values for all of the LC modules configured for the instrument.



Method Editor

Comprehensive Pesticides dMRM\_2.1x150mm\_mod3.m

Properties DA Multisampler Multisampler Pretreatment Binary Pump Column Comp. TQ

Binary Pump (G7112B)

Flow: 0.100 mL/min

Solvents

A: 1 H2O +FA  
2 H2O

B: 1 H2O  
2 ACN

Pressure Limits

Min: 0.00 bar Max: 600.00 bar

Sloptime:  As Injector/No Limit  Off

Posttime:  1.00 min  1.00 min

Advanced

Timetable (empty)

function centric view

Time [min]	A [%]	B [%]	Flow [mL/min]	Max. Pressure Limit [bar]
0.00	50.0	50.0	0.100	600.00

Add Remove Clear All Clear Empty

Cut Copy Paste Shift Times 0.00 min

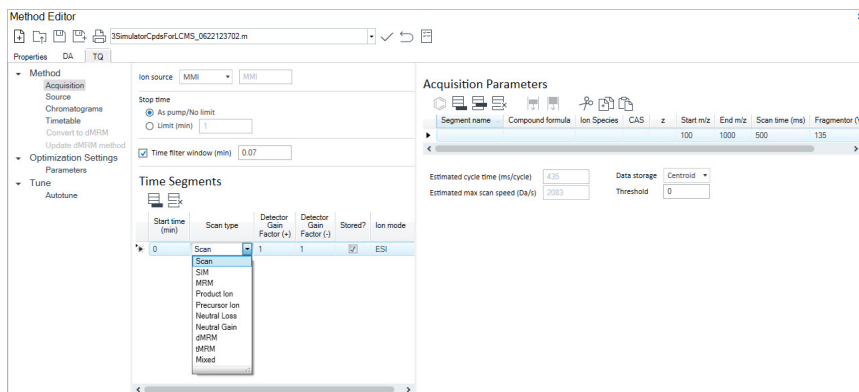
- 3 Enter the TQ parameter values.
  - a On the Acquisition tab, select the Scan type from the list in the **Time Segments** table. The Acquisition Parameters table is cleared when you change the **Scan type**. The parameters available on the right change depending on the **Scan type** for the selected **Time Segment**.

If you are changing the **Scan type** from **MRM** to **Dynamic MRM** or to **Triggered MRM**, you can copy and paste the transitions from the original **Scan segments** table to the Clipboard and then to the new **Scan segments** table. See the *online Help* for more information.

## Basic Operation

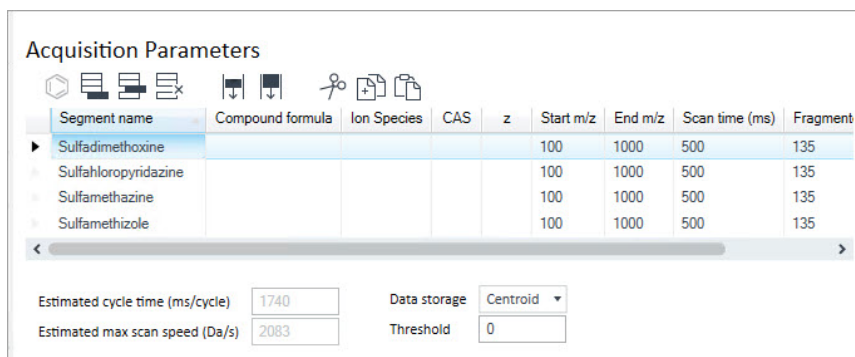
### Setting Up an Acquisition Method

- b Enter any Acquisition values you want to change. Multiple **Scan segments** are allowed.



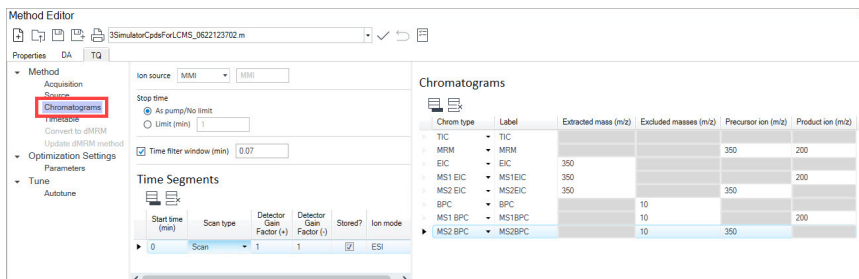
- 4 Set up to change instrument parameters with segments and scans:

- a To add a segment, click **Add Row** or **Insert Row**.
- b Enter the parameters for each Scan segment.



- 5 *Optional.* Set up signals for the Chromatogram plot:

- a Click **Chromatograms**.
- b Select the Chrom Type, and type other plot values.



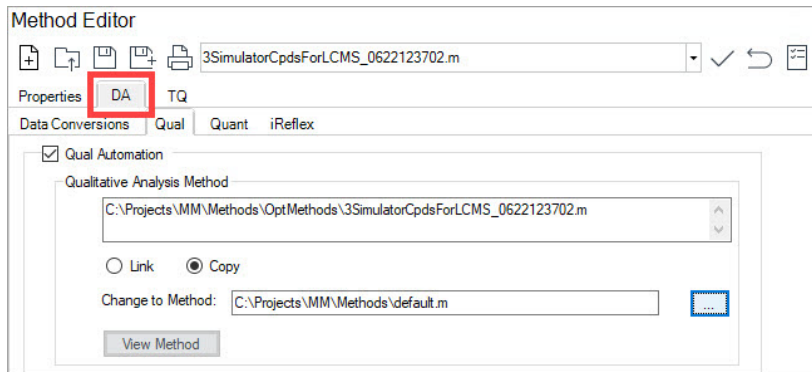
## Basic Operation

### Setting Up an Acquisition Method

#### 6 Set up the data analysis (DA) parameters.

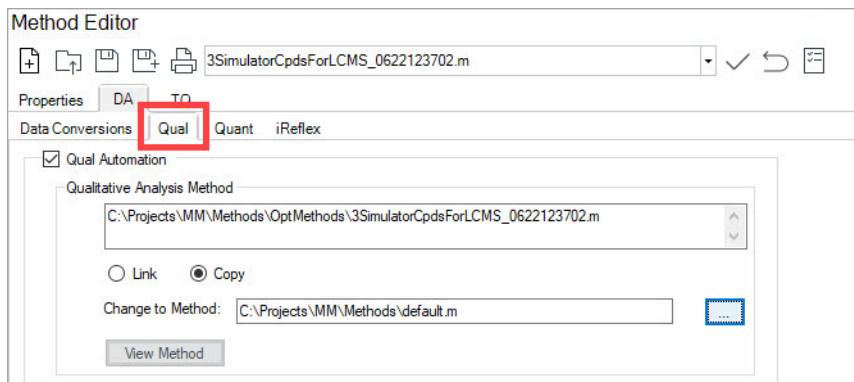
A method can contain Qualitative Analysis parameters, Quantitative Analysis parameters, or a combination. The **Qual Automation** check box is marked on the Qual tab and the **Quant Automation** check box is marked on the Quant tab.

##### a Select the DA tab.

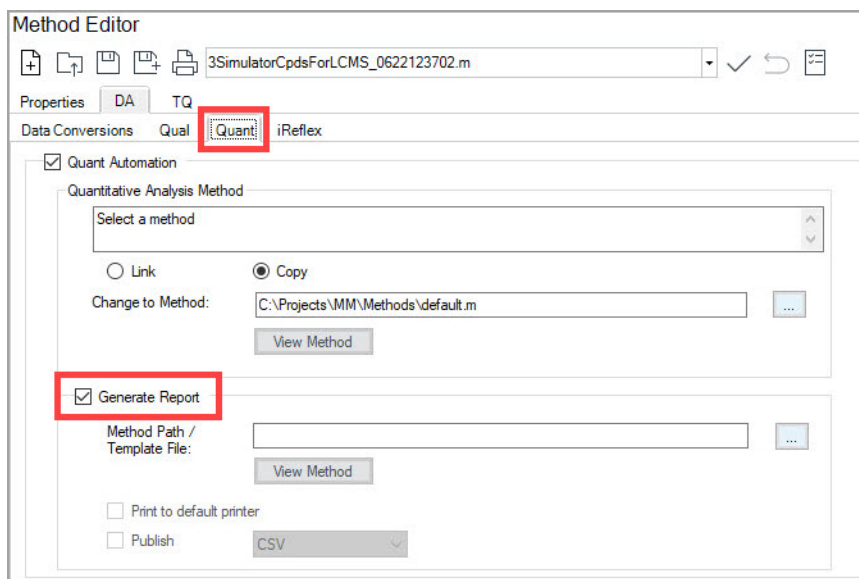


##### b *Optional.* For information on the **Data Conversions** tab, refer to the online Help.

##### c *Optional.* On the **Qual** tab, mark the **Qual Automation** check box. The name of the current Qualitative Analysis method is shown in the box. If you want to change the Qualitative Analysis method that is connected, click **browse** to select a different method. When the Acquisition method is saved, the Qualitative Analysis method that you selected is copied or linked to the Acquisition method.



- d *Optional.* On the **Quant** tab, mark the **Quant Automation** check box. The name of the current Quantitative Analysis method is shown in the list. If you want to change the Quantitative Analysis method that is connected, click **...** **Browse** to select a different method. When the Acquisition method is saved, the Quantitative Analysis method that you selected is copied or linked to the Acquisition method.
- e *Optional.* Mark the **Generate Report** check box on the Quant tab. Then, you select the **Method Path/Template File** to use. If you want to print the report, mark the **Print to default printer** check box. Mark the **Publish** check box to create a CSV file, a TXT file, or a PDF file.



- f *Optional.* For information on the **iReflex** tab, refer to the online Help.
- 7 *Optional.* On the **Properties** tab, set up the properties for this method.
- 8 Save the method.
  - a Click **Save** or **Save As**.
  - b If the Audit Trail Settings in Control Panel are set to prompt for a reason when the method is saved, enter the **Reason for creating a new version of this method**. Refer to the MassHunter Control Panel online Help for more information. Click **OK**.
  - c If necessary, name the method and click **OK**.

# Setting Up and Running a Single Sample

- 1 Click the **Sample Run** window.
- 2 Enter the **Sample Name**, the **Data File Name** and **Path**, and other values.
- 3 *Optional.* Enter the **Additional Information**. If needed, change the value of the parameters in the **Additional Information** list.

Sample Run

Sample

Name:  Position:

Injection Volume:  µL

Comment:

Data File

Auto Increment

Name:

Path:

Additional Information

Parameter Name	Parameter Value
Sample ID	
Override DA Method	
Part of Method to run	Acquisition Only
Equilib Time (min)	0


Method Editor | **Sample Run** | Worklist

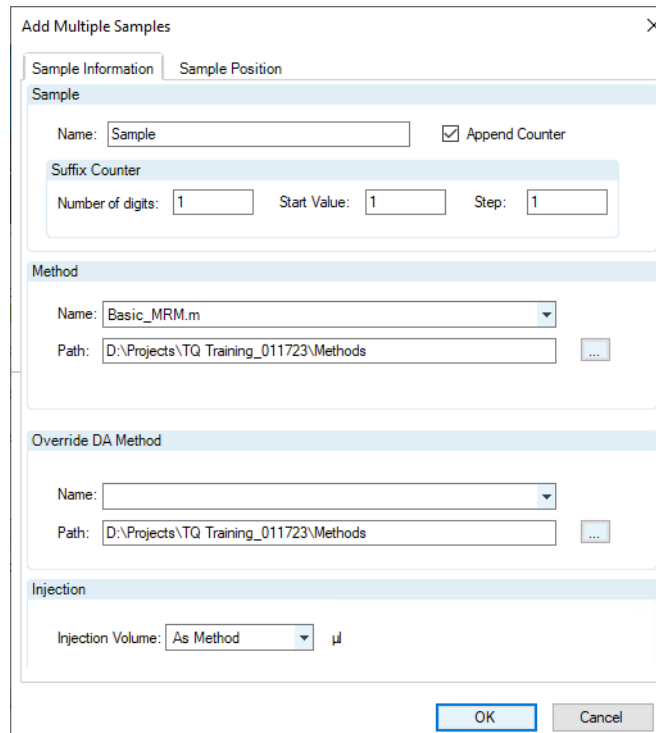
Run a Data Analysis method from this window by selecting **Both Acquisition and DA** or **Acquisition Only** for the **Part of Method to run**. In addition, specify the **Override DA method** to indicate which Acquisition method contains the DA (Data Analysis) method to execute. You always have to do this.

- 4 To start the single sample run, click ► **Run** in the Sample Run window toolbar.



## Setting Up and Running a Worklist

- 1 Click **Worklist** to show the Worklist window.
- 2 Click **Add Multiple Samples** . The **Add Multiple Samples** dialog box opens.
- 3 Enter all the information on the **Sample Information** tab.



The screenshot shows the 'Add Multiple Samples' dialog box with the 'Sample Information' tab selected. The dialog has a title bar with a close button (X) and two tabs: 'Sample Information' and 'Sample Position'. The 'Sample Information' tab contains the following fields:

- Sample** section:
  - Name:
  - Append Counter
- Suffix Counter** section:
  - Number of digits:
  - Start Value:
  - Step:
- Method** section:
  - Name:
  - Path:  ...
- Override DA Method** section:
  - Name:
  - Path:  ...
- Injection** section:
  - Injection Volume:   $\mu$


At the bottom right, there are 'OK' and 'Cancel' buttons.

## Basic Operation

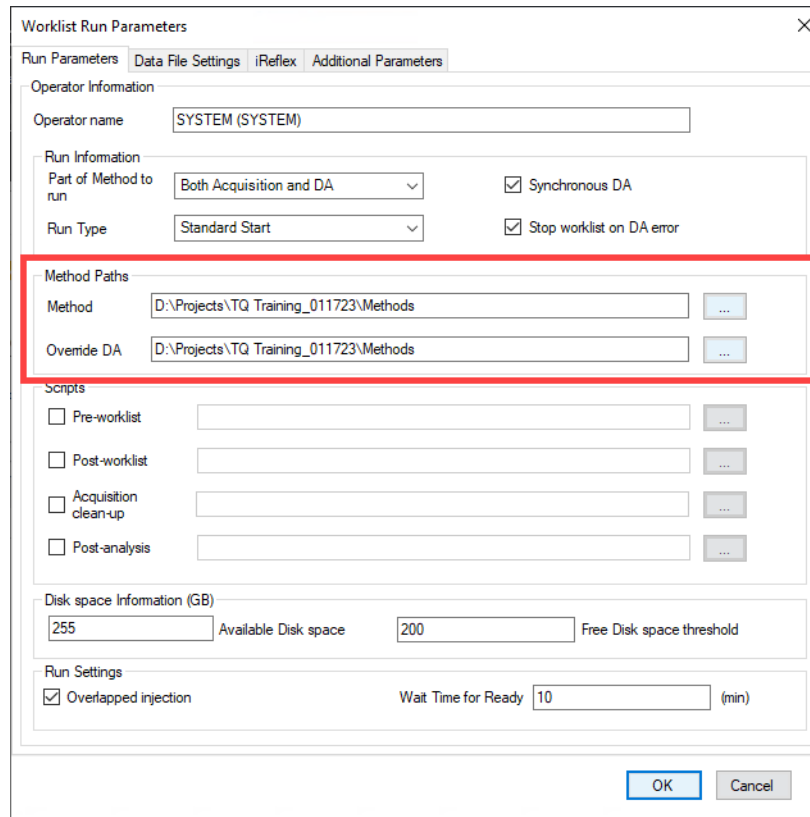
### Setting Up and Running a Worklist

- 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).

The screenshot shows the 'Add Multiple Samples' dialog box with the 'Sample Position' tab selected. The 'Current Configuration' checkbox is unchecked. Under 'Select Well-plate or Vial Tray', a tree view shows 'Well-plate Tray' expanded with 'Plate 1' selected. The 'Plate/Tray Type' dropdown is set to '\*96Agilent\*'. The 'Selection Origin' section has 'Top left' selected. The 'Block Increment' section has 'Row major' selected. The 'Number of samples' field contains '0' and the 'Number of replicates' field contains '1'. Below these settings is a 12x8 grid of wells labeled A-H and 1-12. The grid is highlighted with a red border. At the bottom are 'OK' and 'Cancel' buttons.

- Specify the locations, and click **OK**.
- To set up the worklist run, click **Worklist Run Parameters** .

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



Worklist Run Parameters

Run Parameters | Data File Settings | iReflex | Additional Parameters

Operator Information

Operator name: SYSTEM (SYSTEM)

Run Information

Part of Method to run: Both Acquisition and DA (dropdown)  Synchronous DA

Run Type: Standard Start (dropdown)  Stop worklist on DA error

Method Paths

Method: D:\Projects\TQ Training\_011723\Methods (dropdown) ...

Override DA: D:\Projects\TQ Training\_011723\Methods (dropdown) ...

Scripts

Pre-worklist (dropdown) ...

Post-worklist (dropdown) ...

Acquisition clean-up (dropdown) ...

Post-analysis (dropdown) ...

Disk space Information (GB)

255 Available Disk space 200 Free Disk space threshold

Run Settings

Overlapped injection Wait Time for Ready 10 (min)


OK Cancel

- 8 On the **Data File Settings** tab, enter or select the folders for the data files. Select the **File Naming** options.

The screenshot shows the 'Worklist Run Parameters' dialog box with the 'Data File Settings' tab selected. The dialog has four tabs: 'Run Parameters', 'Data File Settings', 'iReflex', and 'Additional Parameters'. The 'Data File Settings' tab is active and contains the following sections:

- Root Folder:** A text field containing 'D:\Projects\TQ Training\_011723\Data' and a browse button (...).
- Sub-folders:** Three dropdown menus labeled 'Sub-folder 1', 'Sub-folder 2', and 'Sub-folder 3', all currently set to '<Empty>'. Each dropdown has an upward-pointing arrow on its right side.
- File Naming:** Three dropdown menus labeled 'Part 1', 'Part 2', and 'Part 3'. 'Part 1' is set to 'User Text', 'Part 2' is set to 'Counter[0001]', and 'Part 3' is set to '<Empty>'. Below these is a text field containing 'WorklistData'. A checked checkbox labeled 'Use separator for file name parts' is followed by a 'Separator:' label and a text field containing '-'. Each dropdown menu has an upward-pointing arrow on its right side.
- Tree View For File Path and Name:** A tree view showing the file path 'D:\Projects\TQ Training\_011723\Data' and a sub-item 'WorklistData-0001.d'.

At the bottom right of the dialog are 'OK' and 'Cancel' buttons.

- 9 For information on the **iReflex** tab, refer to the online Help.
- 10 *Optional.* On the **Additional Parameters** tab, enter a comment, and click **OK**.
- 11 To start the worklist, click  **Run Worklist**.

#### NOTE

To use an acquisition method that has a different data analysis (DA) method than the method entered in the worklist, show the column called **Override DA Method** in the worklist by using the **Show/Hide/Order Columns** dialog box. In this column, browse for and select the method containing the DA parameters you want to use for the sample. The DA part of this method is used instead of the DA part of the current method.

Or, select this method in the **Add Multiple Samples** dialog box.

## Reviewing Results with Qualitative Analysis

Use the Qualitative Analysis program to do these tasks and more:

- Review results for acquisition method development
- Select the most appropriate precursor and product ions for MRM analyses
- Find compounds
- Identify compounds
- Do molecular feature extraction

Refer to the online Help for the Qualitative Analysis software to learn more.

## Analyzing Data with Quantitative Analysis

Another primary tool for analyzing and reporting instrument results is the Quantitative Analysis program.

Do the exercises in the *Quantitative Analysis Familiarization Guide* to learn how to quantitate the acquired data files:

- Set up a batch and a method to automatically quantitate a set of samples
- Review results by learning how to view and use the Batch-at-a-Glance results screen
- Identify and use outliers to change the method and requantitate the data using a better calibration curve fit or other more appropriate settings

Refer to the online Help for the Quantitative Analysis program to learn how to do more operations to analyze your data.

# 5

## Maintenance

Routine Tasks 64

Calibrant Delivery System and Divert Valve 64

Sources 65

Nebulizer 68

Periodic Tasks 69

Calibrant Delivery System and Divert Valve 69

Sources 70

Nebulizer 75

Corona Needle 80

Ion Injector 82

Rough Pump 85

Disinfection 90

When performing maintenance, it is recommended that you track procedures and details in a maintenance logbook.

## Routine Tasks

### **Calibrant Delivery System and Divert Valve**

The following procedures describe routine maintenance tasks for the calibrant delivery system (CDS) and divert valves.

#### **Checking calibrant levels**

Perform this procedure monthly or weekly if you tune the instrument frequently.

Enough tuning mix must be present to immerse the end of the intake tube.

If the tuning mix level is within a few millimeters of the end of the intake tube, refill the calibrant bottle.



## Sources

The following procedures describe routine maintenance for the components that make up the source.

### Cleaning the spray chamber - daily

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

#### NOTE

Use the weekly cleaning procedure if symptoms of contamination persist or if the spray shield or capillary cap show significant discoloration that cannot be removed by the regular, daily cleaning.

Tools needed:

- Clean powder-free nitrile gloves
- Clean lint-free cloth
- LC/MS-grade isopropanol
- LC/MS-grade water
- Clean wash bottle

To perform daily cleaning of the spray chamber,

- 1 Put on clean powder-free nitrile gloves.
- 2 Lower the drying gas temperature to the minimum level.
- 3 For APCI/multimode, lower the vaporizer heater temperature to the minimum level.

#### WARNING

**Make sure the source is cooled before you clean the spray shield.**

- 4 In a clean wash bottle, prepare a 50:50 mix of LC/MS-grade isopropanol and LC/MS-grade water.
- 5 Open the spray chamber.
- 6 Dampen a clean lint-free cloth with a mixture of LC/MS-grade isopropanol and LC/MS-grade water.
- 7 Wipe the spray shield and the area around the spray shield.
- 8 Close the spray chamber.

**Cleaning the spray chamber - weekly**

Perform this maintenance weekly or whenever symptoms indicate that contamination exists in the spray chamber and normal daily cleaning does not correct the problem.

Tools needed:

- Clean powder-free nitrile gloves
- Clean lint-free cloth
- LC/MS-grade isopropanol
- LC/MS-grade water
- Clean wash bottle
- 4000 grit abrasive paper

To perform weekly cleaning of the spray chamber,

- 1** Put on clean powder-free nitrile gloves.
- 2** Disconnect the nebulizer sample line and sample nebulizer gas tubing from the nebulizer.
- 3** Disconnect all cables and tubings from the instrument.
- 4** Open the latch of spray chamber.
- 5** Open the spray chamber.
- 6** Put the source under a fume hood.
- 7** Dampen a clean lint-free cloth with a mixture of LC/MS-grade isopropanol and LC/MS-grade water. For Multimode, be careful not to touch the thermocouple probe.
- 8** Wipe the interior of the spray chamber with a clean lint-free cloth.
- 9** Make sure you wear clean gloves. Use a tech wipe to hold the parts as you remove them. They will be warm.
- 10** Clean the spray shield, capillary cap, end plates, and contact ring under fume hood. Carefully remove the canted coil spring from the capillary cap before cleaning.
- 11** Use 4000 grit abrasive paper to clean contaminants from the parts.
- 12** Put the spray shield, capillary cap, end plates, and contact ring in the beaker.
- 13** Add a 50:50 mix of LC/MS-grade isopropanol and LC/MS-grade water in a beaker.
- 14** Put the beaker in sonication for 5 minutes.
- 15** Drain the solvent and remove the parts from the beaker.

- 16 Dry the parts by using nitrogen gas.
- 17 Carefully insert the canted coil back into the capillary cap.
- 18 Reinstall the contact ring, end plate, capillary cap, and spray shield.
- 19 Reinstall the spray chamber on the instrument.
- 20 Close the spray chamber.
- 21 Close the latch of spray chamber.
- 22 Connect all cables and tubings to the instrument.

**CAUTION**

**Do not bend or blunt the tip of the needle, as it will decrease system performance. Sharpening the needle is not needed.**

---

- 23 Pull and twist the 4000 grit abrasive paper along the needle and off the tip of the needle.

**CAUTION**

**Do not hit the tip of the corona needle. The tip can bend, which can decrease system performance.**

---

- 24 Starting at the base of the needle, wipe the needle with a clean lint-free cloth. The cloth can be dry or dampened with LC/MS-grade isopropanol.

**CAUTION**

**Do not hit the tip of the corona needle. The tip can bend, which can decrease system performance.**

---

- 25 Connect the nebulizer sample line and sample nebulizer gas tubing to the nebulizer.

## Nebulizer

The following procedures describe routine maintenance for the nebulizer.

### Flushing the nebulizer

Check the nebulizer weekly. If needed, flush traces of samples and buffers out of the tubing, valves, and nebulizer.

Parts needed:

- LC/MS-grade methanol

To flush the nebulizer,

- 1** Install a mix of 30% to 70% methanol in water.  
On a monthly basis, prepare enough mix to flush overnight.
- 2** On a daily basis, flush for 3 minutes.  
On a monthly basis, flush overnight.

## Periodic Tasks

### Calibrant Delivery System and Divert Valve

The following procedures describe periodic maintenance tasks for the calibrant delivery system (CDS) and divert valves.

#### Filling a calibrant bottle

Perform this procedure as needed.

Parts required:

- APCI calibrant
- APCI-L tuning mix
- ESI-L tuning mix
- MMI-L tuning mix
- Reference mix

- 1 Rinse the calibrant bottle before refilling.
- 2 Refill the calibrant bottle with the appropriate tuning mix.
- 3 Put the intake tube into the calibrant bottle as you lift the calibrant bottle into position.

#### CAUTION

**Do not touch the intake tube unless you are wearing clean gloves.**

- 4 Attach the calibrant bottle onto the fixed bottle cap. Turn the calibrant bottle counter clockwise to tighten.

#### CAUTION

**The bottle only needs to be snug. Do not over-tighten the bottle. A leaky bottle cap can prevent the bottle from pressurizing and can lead to the evaporation of the contents.**

#### CAUTION

**The tuning mixes are not interchangeable. Failure to change the tuning mix when the source or inlet assembly is changed can result in miscalibration of the instrument and errors in mass assignments.**

## Sources

The following procedures describe periodic maintenance for the components that make up the source. See **Table 2** for tuning mixes.

### Changing to a source that uses a different tuning mix

Complete these steps every time you change a source that uses a different tuning mix.

- 1 Right-click the TQ device pane and select **Calibrant**.
- 2 Run the calibrant for 5 minutes.
- 3 Perform a checktune. See **“Calibrating the LC/TQ (Checktune)”** on page 29.
- 4 If the checktune results are not acceptable, run autotune on a supported tuning source. See **“Calibrating the LC/TQ (Autotune)”** on page 30.
- 5 When the instrument is ready, load or create the method to use with the source.
- 6 Change the method settings for the new source.

### Removing an Electrospray or Agilent Jet Stream source

Before you begin:

- Do a complete autotune (see **“Calibrating the LC/TQ (Autotune)”** on page 30). If an autotune is not available, do a checktune (see **“Calibrating the LC/TQ (Checktune)”** on page 29).
- If the tune report does not show good results, tune the instrument. If needed, change the source to one that supports autotune.
- Put the system in Standby mode.

#### NOTE

In the Tune view or context, turn down the gas temperature as low as possible. Leave the gas flow on for 30 minutes to cool the source.

- 1 For Electrospray (ESI), disconnect the nebulizer sample line.
- 2 For the dual ESI or dual AJS source on TOF and Q-TOF systems only, remove the second nebulizer. Disconnect the nebulizer sample line and sample nebulizer gas tubing.
- 3 Disconnect the heating cable.
- 4 Disconnect the sheath gas line at the connector and unplug it.

- 5 Disconnect the AJS HV cable delivering AJS nozzle voltage.
- 6 Disconnect the Multimode cable delivering AJS nozzle voltage.
- 7 Open the latch of spray chamber.
- 8 Open the spray chamber.
- 9 Remove the spray chamber by lifting it along the axis of its hinges.

**WARNING**

**The spray shield can be hot. Be careful not to burn yourself when you remove the spray shield.**

- 10 If you are changing to a different source type, remove the spray shield.

**Removing an APCI source**

- 1 Put the system in Standby mode.

**NOTE**

In the Tune view or context, turn down the gas temperature as low as possible. Leave the gas flow on for 30 minutes to cool the source.

- 2 Disconnect the nebulizer sample line.
- 3 Disconnect the sample nebulizer gas tubing.
- 4 Disconnect the heater cable.
- 5 Disconnect the corona needle voltage cable.

**Removing a Mutlimode source**

- 1 Put the system in Standby mode.

**NOTE**

In the Tune view or context, turn down the gas temperature as low as possible. Leave the gas flow on for 30 minutes to cool the source.

- 2 Disconnect the nebulizer sample line.
- 3 Disconnect the heater cable.
- 4 Disconnect the charging electrode cable.
- 5 Open the latch of spray chamber.

**WARNING**

The spray chamber operates at very high temperature. Do not continue until the spray chamber is cool.

---

- 6 Remove the spray chamber by lifting it along the axis of its hinges.

**WARNING**

The spray shield can be hot. Be careful not to burn yourself when you remove the spray shield.

---

- 7 If you are changing to a different source type, remove the spray shield.

**Installing an Electrospray or Agilent Jet Stream source**

- 1 Check that the appropriate capillary cap is installed.
- 2 Install the appropriate spray shield.
- 3 For Electrospray (ESI), do the following:
  - a Make sure the small hole is at the top ("12 o'clock position").
  - b If needed, use a T10 Torx screwdriver to loosen the two screws in the end plate.
  - c Rotate the ESI spray shield clockwise until the hole is in the correct position.
  - d Gently tighten the Torx screws again.
- 4 Install the spray chamber.
- 5 Close the spray chamber.
- 6 Connect the MultiMode cable delivering AJS nozzle voltage.
- 7 Connect the heater cable.
- 8 Connect the sample nebulizer gas tubing.
- 9 Connect the nebulizer sample line.

**CAUTION**

Do not hit the tip of nebulizer needle. The needle is not protected. If you damage the needle, system performance can be affected.

---

- 10 Connect the reference nebulizer sample line and sample nebulizer gas tubing.  
In the Acquisition program,



- 1 Turn on the LC/MS instrument.
- 2 While you wait for the set points to equilibrate, prepare and install calibrant for the new ion source.
  - a Rinse a clean calibrant bottle with LC/MS-grade acetonitrile.
  - b Pour the ESI Calibrant into the calibrant bottle.
  - c Install the calibrant bottle into the calibrant delivery system.

### Installing an APCI source

- 1 Install the appropriate spray shield.
- 2 Make sure the small hole is at the top ("12 o'clock position").
- 3 If needed, use a T10 Torx screwdriver to loosen the two screws in the end plate.
- 4 Rotate the ESI spray shield clockwise until the hole is in the correct position.
- 5 Gently tighten the Torx screws again.
- 6 Install the spray chamber.
- 7 Close the spray chamber.
- 8 Close the spray chamber latch.
- 9 Connect the heater cable.
- 10 Connect the corona needle voltage cable.
- 11 Connect the sample nebulizer gas tubing.
- 12 Connect the nebulizer sample line.

In the Acquisition program,

- 1 Turn on the LC/MS instrument.
- 2 While you wait for the set points to equilibrate, prepare and install calibrant for the new source.
  - a Rinse a clean calibrant bottle with LC/MS-grade acetonitrile.
  - b Pour the APCI Calibrant into the calibrant bottle.
  - c Install the calibrant bottle into the calibrant delivery system.

### Installing a Mutlimode source

- 1 Install the appropriate spray shield.
- 2 Make sure the dowels are at 6 and 9 o'clock positions.
- 3 Install the spray chamber.

- 4 Close the spray chamber.
- 5 Close the latch of spray chamber.
- 6 Connect the corona needle voltage cable.
- 7 Connect the charging electrode cable.
- 8 Connect the heater cable.
- 9 Connect the sample nebulizer gas tubing.
- 10 Connect the nebulizer sample line.

In the Acquisition program,

- 1 Turn on the LC/MS instrument.
- 2 While you wait for the set points to equilibrate, prepare and install calibrant for the new source.
  - a Rinse a clean calibrant bottle with LC/MS-grade acetonitrile.
  - b Pour the MMI Calibrant into the calibrant bottle.
  - c Install the calibrant bottle into the calibrant delivery system.

## Nebulizer

The following procedures describe periodic maintenance for the nebulizer.

### Removing the nebulizer

Perform this procedure when you need to access the nebulizer for maintenance.

Tools needed:

- Clean powder-free nitrile gloves

To remove the nebulizer,

- 1 Put on clean powder-free nitrile gloves.
- 2 In the Acquisition program, shut off the LC solvent flow and nebulizing gas.
- 3 Remove the left side cover.
- 4 Disconnect the nebulizer sample line from the nebulizer.
- 5 Disconnect the sample nebulizer gas tubing from the nebulizer.

### WARNING

**The tip of the nebulizer can be very hot. Do not touch the nebulizer until it is cool.**

- 6 Turn the nebulizer counterclockwise until it disengages from the retaining screws.
- 7 Gently lift the nebulizer out of the spray chamber.

### Inspecting the nebulizer

Perform this procedure if you have the G1960-67470 Nebulizer Adjustment Kit. Perform this project if there is increased LC back pressure, off-axis spraying, dripping from the nebulizer, or when the reference nebulizer needle is not spraying.

Tools needed:

- Nebulizer Adjustment Fixture
- Clean powder-free nitrile gloves
- T6 Torx driver
- 1/4-inch x 5/16-inch open-end wrench (Quantity: 2)

Parts needed:

- Replacement needle kit

To inspect the nebulizer,

- 1 Put on clean powder-free nitrile gloves.
- 2 Determine your nebulizer type.
- 3 Carefully insert the tip of the nebulizer into the end of the Nebulizer Adjustment Fixture.
- 4 Slide the heads of the two spring loaded socket head cap screws through the mounting holes on the body of the nebulizer.
- 5 Turn the nebulizer body 1/8 turn counterclockwise.

**NOTE**

If the nebulizer is too tight or too loose, loosen or tighten the nut on the captured screws to adjust the tension on the spring. Turn clockwise to tighten, counterclockwise to loosen.

- 6 Gently push the 40 x Magnifier into the end horizontal position of the Nebulizer Adjustment Fixture. Align the tapered tip of the magnifier to the tip of the nebulizer needle.
- 7 Push the LED Flashlight into the Nebulizer Adjustment Fixture. No tool is needed.
- 8 Turn on the LED Flashlight. The on/off switch is at the end of the flashlight.

**CAUTION**

**Do not hit or touch the end of the nebulizer or you can damage the needle.**

- 9 Gently push the LED Flashlight into the end horizontal position of the Nebulizer Adjustment Fixture.
- 10 Push the 40 x Magnifier into the Nebulizer Adjustment Fixture. No tool is needed. Align the tapered tip of the magnifier to the tip of the nebulizer needle.
- 11 Turn on the LED Flashlight. The on/off switch is at the end of the flashlight.

**CAUTION**

**Do not hit or touch the end of the nebulizer or you can damage the needle.**

- 12** Gently slide and rotate the 40 x Magnifier into or out of the Nebulizer Adjustment Fixture to focus the magnifier on the end of the nebulizer nozzle and needle. For the APCI nebulizer, if needed, remove the pocket clip to focus properly.

Refer to **Table 3** for correct needle position.

**Table 3** Correct needle position

Nebulizer	Distance
Microflow Nebulizer Assembly	0.003 inch
ES Tested Nebulizer	0.000 (flush)
Tested ESI Nebulizer Assembly (older style)	0.003 inch
Tested APCI Nebulizer Assembly	0.000 (flush)

### Adjusting or replacing the nebulizer needle

Perform this procedure if you have the G1960-67470 Nebulizer Adjustment Kit. Perform this project if there is increased LC back pressure, off-axis spraying, dripping from the nebulizer, or when the reference nebulizer needle is not spraying.

Tools needed:

- Nebulizer Adjustment Fixture
- Clean powder-free nitrile gloves
- T6 Torx driver
- 1/4-inch x 5/16-inch open-end wrench (Quantity: 2)

Parts needed:

- Replacement needle kit

To adjust or replace the nebulizer,

- 1 Put on clean powder-free nitrile gloves.
- 2 Determine the nebulizer holder type.
- 3 Install the nebulizer in the Nebulizer Adjustment Fixture.
- 4 Loosen the locknut next to the zero-dead-volume (ZDV) union.
- 5 Remove the union from the nebulizer needle holder.
- 6 Remove the union lock nut from the nebulizer.

- 7 Loosen the fastener (locknut, 1.5-mm hex screw or T6 Torx screw) that secures the needle holder in place.
- 8 Loosen the needle holder and pull it out of the nebulizer body. Dispose of the used needle properly.
- 9 For more thorough cleaning, sonicate the nebulizer body. Pour isopropyl alcohol or methanol through inside of the shaft.
- 10 Thread the new needle holder into the nebulizer body.
- 11 Carefully remove the new needle from the shipping tube and adhesive strip. Identify the sharp end of the needle.
- 12 Carefully slide the sharp end of the needle through the ferrule until 1 cm of the needle remains at the ferrule.

**CAUTION**

**Do not bend the needle as you slide it into the ferrule. The needle is fragile. A bent needle results in poor alignment with the nebulizer body.**

- 13 Push the needle into the ferrule until the needle is slightly visible flush with the ferrule.
- 14 Reinstall the locknut and the union. Hand-tighten the union.
- 15 Hold the needle holder steady with a 3-mm open-end wrench or 1/4-inch x 5/16-inch open-end wrench, depending on your nebulizer holder style.
- 16 Tighten the union one-half to three quarters of a turn, or until an audible squeak is heard.  
The ferrule is now compressed.
- 17 Pull carefully on the needle to ensure the needle is held firmly in place.
- 18 Insert the needle into the nebulizer shaft.
- 19 Adjust the electrospray needle position before reinstalling the nebulizer in the spray chamber.
- 20 Tighten the fastener (T6 Torx screw) that secures the needle holder in place.

### Installing the nebulizer

- 1 Insert the nebulizer part way into the spray chamber.

**CAUTION**

**Do not hit the tip of the needle as you insert the nebulizer. The tip of the needle is easily damaged.**

- 2 Connect the nebulizer sample line to the nebulizer.
- 3 Turn the nebulizer clockwise and lock it in place.
- 4 Connect the sample nebulizer gas tubing to the nebulizer.

**CAUTION**

**Do not over-tighten the LC fitting. An over-tightened fitting can crush the tubing, which creates a restriction.**

### Fixing Spray Stability EMF

If spray stability EMF is active, you should inspect, adjust, or flush your nebulizer to ensure a stable spray from the CDS to the nebulizer. An unstable spray may lead to issues with tuning and sample analysis.

- 1 Inspect ([page 75](#)), adjust ([page 77](#)), or flush ([page 68](#)) your nebulizer. If necessary, replace the nebulizer needle or nebulizer ([page 77](#)).
- 2 Follow the procedure **“Calibrating the LC/TQ (Checktune)”** on page 29 or **“Calibrating the LC/TQ (Autotune)”** on page 30. The spray stability is reevaluated as part of a tune. If it is within the acceptable range, the EMF warning should be removed following completion of the tune.

## Corona Needle

The following procedures describe periodic maintenance for the corona needle.

### Cleaning the corona needle

Perform this procedure when you observe decreased sensitivity, decreased signal stability, or increased corona voltage during APCI operation.

Tools needed:

- Clean powder-free nitrile gloves
- Clean lint-free cloth
- LC/MS-grade isopropanol
- 4000 grid abrasive paper

To clean the corona needle,

1 Put on clean powder-free nitrile gloves.

### WARNING

**The needle and related parts can be very hot. Do not touch the needle or related parts until they are cool.**

2 Pull the corona needle assembly out of the spray chamber.

### CAUTION

**Do not bend or blunt the tip of the needle, it will decrease system performance. Sharpening the needle is not needed.**

3 Fold a piece of 4000 grit abrasive paper over the base of the needle.

4 Pull and twist the 4000 grit abrasive paper along the needle and off the tip of the needle.

### CAUTION

**Do not hit tip of the corona needle. The tip can bend, which can decrease system performance.**

5 Starting at the base of the needle, wipe the needle with a clean lint-free cloth. The cloth can be dry or dampened with LC/MS-grade isopropanol.

6 Reinstall the corona needle assembly in the spray chamber.



**Replacing the corona needle**

Perform this procedure when symptoms indicate poor corona needle performance and cleaning the needle does not restore performance.

**CAUTION**

**Do not hit the tip of the corona needle. The tip can bend, which can decrease system performance.**

Tools needed:

- Clean powder-free nitrile gloves
- Corona needle

To replace the corona needle,

- 1 Put on clean powder-free nitrile gloves.
- 2 Pull the corona needle assembly out of the spray chamber.
- 3 Remove the needle collar from the corona needle shaft.
- 4 Remove the old corona needle from the collar.
- 5 Install a new corona needle, with its integral ferrule, in the collar.
- 6 Turn the collar onto the needle holder and tighten by hand.

**CAUTION**

**Do not hit the tip of the corona needle. The tip can bend, which can decrease system performance.**

- 7 Reinstall the corona needle assembly in the spray chamber.

## Ion Injector

The following procedures describe periodic maintenance for the ion injector. These procedures apply to both single and multi-bore ion injectors.

### Removing the ion injector

Perform this procedure when you need to clean or replace the ion injector.

Tools needed:

- Clean powder-free nitrile gloves
- Capillary Puller Tool

#### WARNING

**The spray chamber operates at very high temperatures. Do not continue until the spray chamber is cool.**

To remove the ion injector,

- 1 Put on clean powder-free nitrile gloves.
- 2 Remove the ion injector cap from the end of the ion injector.
- 3 Screw the capillary puller tool fully into the spray shield mount.
- 4 Push the collet assembly over the ion injector and then tighten by holding the puller handle. Turn the collet clamping knob clockwise.
- 5 Pull on the knob until the collet assembly stops moving.
- 6 Visually check that the ion injector has been pulled out by the collet by looking through the slots on the tool.
- 7 Rotate the knob assembly counterclockwise slightly, and pull out the assembly to unlock the collet assembly from the puller.
- 8 Pull the tool out.

#### CAUTION

**Carefully pull the ion injector out along its long axis. The ion injector is glass or of similar material, and you can break it by putting vertical or horizontal pressure on it.**

- 9 Loosen the collet clamping knob slightly to loosen the ion injector from the capillary puller tool.
- 10 Pull out the ion injector from the collet.
- 11 Unscrew the tool from the shield mount.

**Cleaning the ion injector**

Perform this procedure when you observe decreased sensitivity and decreased signal stability.

Tools needed:

- Clean powder-free nitrile gloves
- LC/MS-grade isopropanol
- Alconox powdered precision cleaner
- Deionized (18 M $\Omega$ /cm) water
- 100-mL polypropylene graduated cylinder
- 10-mL polypropylene graduated cylinder
- Two 1-mL pipette tip

To clean the ion injector,

- 1 Put on clean powder-free nitrile gloves.
- 2 Dissolve 1 gram of Alconox Powdered Precision Cleaner in 100 mL of deionized (18 M $\Omega$ /cm) water. This concentration is the recommended concentration for both manual or ultrasonic cleaning.
- 3 If you use a glass graduated cylinder, cover the end of the ion injector with a 1-mL pipette tip to protect the metalized plating.
- 4 Trim the pipette tip to approximately 4 cm so that the ion injector can be immersed in the cleaning solution.
- 5 Place the ion injector upright in a graduated cylinder and fill with Alconox Powdered Precision Cleaner solution.
- 6 Put the Ultrasonic cleaner under a fume hood. Sonication should be done under a fume hood.
- 7 Sonicate the graduated cylinder that contains the ion injector in an ultrasonic cleaner for 10 to 15 minutes. Do not clean longer than 15 minutes.

**NOTE**

Rinse the ion injector and graduated cylinder several times with deionized water.

- 8 Fill the graduated cylinder with deionized water and sonicate for 10 to 15 minutes.
- 9 Remove the ion injector from the graduated cylinder and remove the pipette tips (if any were used).
- 10 Blow out excess water from the ion injector bore with nitrogen.

**Installing the ion injector**

Perform this procedure after cleaning the ion injector or when installing a new ion injector.

Tools needed:

- Clean powder-free nitrile gloves
- LC/MS-grade isopropanol

To install the ion injector,

- 1 Put on clean powder-free nitrile gloves.
- 2 Lubricate the ion injector entrance end with LC/MS-grade isopropanol.
- 3 Carefully insert the ion injector straight into the desolvation assembly.

**CAUTION**

**Putting vertical and horizontal pressure on the ion injector can break it. Spray ion injector with isopropanol alcohol to lubricate the ion injector for easier removal.**

- 4 Slide the ion injector straight into the desolvation assembly.

The ion injector must be aligned correctly so that its end fits into a fixed ion injector cap inside desolvation assembly.

- 5 When 2 to 3 cm of the ion injector remains extended from the desolvation assembly, the ion injector rests against the rear contact spring, which slightly restricts further insertion of the ion injector. Slightly increase the insertion force to push the ion injector through the rear contact spring to fully insert the injector.
- 6 Continue to apply pressure until approximately 1 cm remains extended from the desolvation assembly.

**CAUTION**

**Do not twist or turn the ion injector cap when you install it or you can damage the metal plating.**

- 7 Install the contact ring and end plate.
- 8 Install the two T10 Torx screws to hold the end plate.
- 9 Lubricate the ion injector tip with LC/MS-grade isopropanol.
- 10 Install the ion injector cap over the outer end of the ion injector.
- 11 Install the spray shield.
- 12 Close the spray chamber.

## Rough Pump

The following procedures describes periodic maintenance of the vacuum system of the instrument.

### Checking the rough pump oil level

- 1 Check that the oil level is between the marks for Max and Min.
- 2 Check that the pump oil is clear and the color is lighter than amber.
- 3 If the pump oil is dark or full of suspended particles, replace it.

### Adding rough pump oil

Perform this procedure when the pump oil level is low.

Tools needed:

- Clean chemical-resistant gloves
- Safety glasses
- Funnel
- 10 mm hex key
- Paper towel or shop rag

Parts needed

- AFV 60 gold oil

---

**WARNING**

**Wear chemical-resistant gloves and safety glasses (goggles) for your safety.**

---

**WARNING**

**Do not touch the fill cap or pump until they are cool. The fill cap and pump can be dangerously hot.**

---

**WARNING**

**Never add or replace the rough pump oil while the pump is on. Hot oil can splash out and cause harm.**

---

**CAUTION**

**Use only the rough pump oil appropriate for your pump (AFV 60 gold oil for Agilent MS40+). The use of any other oils can substantially reduce pump life and will invalidate the pump warranty.**

To add rough pump oil,

- 1 Put on clean chemical-resistant gloves.
- 2 Place the instrument into Standby mode
- 3 Vent the instrument.
- 4 Allow the rough pump to turn off automatically. This should take about 15 minutes.
- 5 Unplug all power cords from LC/MS instrument.
- 6 Unplug the rough pump.
- 7 Remove the fill cap on the rough pump.
- 8 Using a funnel, add new AFV 60 gold oil until the oil level matches the maximum mark.

**CAUTION**

**Do not fill the oil over the maximum mark beside the oil level window.**

- 9 Reinstall the fill cap.
- 10 Wipe off all excess oil around and underneath the pump.

**Replacing the filter**

Perform this procedure yearly.

Tools needed:

- Pair of pliers

Parts needed:

- Filter
- Filter cap O-ring
- Small cap O-ring

To replace the filter,

- 1 Place the instrument into 'Standby' then 'Vent'.
- 2 Allow the rough pump to turn off automatically. This should take about 15 minutes.
- 3 Unplug all power cords from LC/MS instrument.
- 4 Unplug the rough pump
- 5 Disconnect the exhaust tube.
- 6 Unscrew the filter cap.

**CAUTION**

**The filter cap is spring loaded.**

- 7 Remove the O-ring from the filter cap.
- 8 Remove the old filter.
- 9 Remove the O-ring from the filter.
- 10 Install the new small O-ring on the new filter and install the filter into the pump in the proper orientation.
- 11 The distance of the filter should not be < 2 cm from the outside cap thread.
- 12 Make sure the filter is > 2 cm from the outside cap threads. When properly installed:
  - The filter cap is flush to the pump.
  - The back end of the filter is seated inside the filter mount.
- 13 Install the new large O-ring and filter cap.
- 14 Connect the exhaust tube.
- 15 Plug in the rough pump and all power cords from the LC/MS instrument.

**Replacing the rough pump oil**

Perform this procedure yearly.

Tools needed:

- Clean chemical-resistant gloves
- Safety glasses
- Funnel
- 10 mm hex key
- Paper towel or shop rag
- Small plastic tub

Parts needed:

- AFV 60 gold oil

**WARNING**

**Wear chemical-resistant gloves and safety glasses (goggles) for your safety.**

**WARNING**

**Do not touch the fill cap or pump until they are cool. The fill cap and pump can be dangerously hot.**

**WARNING**

**Never add or replace the rough pump oil while the pump is on. Hot oil can splash out and cause harm.**

**CAUTION**

**Use only the rough pump oil appropriate for your pump (AFV 60 gold oil for Agilent MS40+). The use of any other oils can substantially reduce pump life and will invalidate the pump warranty.**

To replace the rough pump oil,

- 1 Put on clean chemical-resistant gloves.
- 2 Place the instrument into 'Standby' then 'Vent'.
- 3 Allow the rough pump to turn off automatically. This should take about 15 minutes.
- 4 Unplug all power cords from the LC/MS instrument.
- 5 Unplug the rough pump.
- 6 Remove first the fill cap and then open the drain plug.



**CAUTION**

The rough pump is very heavy. Use care if you move or tip the pump.

- 7 Allow the oil to drain completely.
- 8 Reinstall the drain plug.
- 9 Using a funnel, add new AFV 60 gold oil until the oil level matches the maximum mark.

**CAUTION**

Do not fill the oil over the maximum mark beside the oil level window.

- 10 Reinstall the fill cap.

**Replacing fuses or resetting circuit breakers**

Perform this procedure as needed for maintenance.

Tools needed:

- Long flat-blade screwdriver
- 8 Amp fuse
- 12.5 Amp fuse

**WARNING**

Never replace a fuse while the instrument is plugged into the power outlet. You run the risk of an electric shock.

**CAUTION**

If the instrument is plugged in and turned on when you change a fuse, you will damage the instrument.

To replace fuses or reset circuit breakers,

- 1 Unplug the instrument power cord or cords from the power outlet.
- 2 If your LC/MS instrument includes circuit breakers and you need to reset them, press in the circuit breaker buttons.
- 3 Using a flat-blade screw driver, remove the fuse holder of the blown fuse.
- 4 Replace with the appropriate fuse.
- 5 Reinstall the fuse holder.
- 6 Plug in the instrument.

## Disinfection

To prevent the spread of respiratory illness, disinfect your instrument as needed.

Equipment needed:

- Clean powder-free nitrile gloves
- Clean lint-free cloth
- Gown
- Isopropanol
- Mask

To disinfect your instrument,

- 1** Put on personal protective gear (gloves, mask, gown).
- 2** Make sure your instrument is cooled to room temperature.
- 3** Dampen a clean lint-free cloth with the 70:30 isopropyl alcohol:water mix.
- 4** Gently wipe external surfaces to be cleaned using a moistened cloth. Do not allow any liquid to drip into the instrument.
- 5** Wipe the outer surface of cables and their connectors, but do not touch the electrical connections (such as pins and plugs inside the cable or connection port).
- 6** Use one cloth per instrument to prevent cross contamination.
- 7** Dispose of the cloth appropriately.
- 8** Allow all surfaces to completely air-dry. No moisture should be present on the instrument surfaces.
- 9** Discard the gloves and other personal protective equipment or clean them in an approved process, and then wash your hands.

[www.agilent.com](http://www.agilent.com)

©Agilent Technologies, Inc. 2023  
April 2023

DocNo D0025685 Rev. A.01  
DE22440912

