



Agilent MassHunter Workstation

Qualitative and Quantitative Analysis Installation Guide

Notices

Document Identification

D0130987 Rev. A.00

May 2025

Copyright

© Agilent Technologies, Inc. 2025

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 the 13.0 revision or higher of Agilent MassHunter Quantitative and Qualitative Analysis, and compatible with MassHunter Quantitative and Qualitative Analysis, until superseded.

Software Manufacturing



Manufactured for Agilent
Technologies
5301 Stevens Creek Blvd
Santa Clara, CA 95051

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.

In This Guide

Use this guide to install or upgrade MassHunter Qualitative Analysis and Quantitative Analysis.

This guide applies to the installation of:

- MassHunter Qualitative Analysis 13.0 and later
- MassHunter Quantitative Analysis 12.1 and later

Table of Contents

Step 1: Prepare for Installation	6
Step 2: Install the Qualitative Analysis Program	7
Step 3: Install or Upgrade the Quantitative Analysis Program	13
Step 4: Install Quantitative Analysis Supplemental Items	20
Step 5: Configure Excel for MassHunter	23



Installation

This document describes the required steps to install MassHunter Quantitative and Qualitative Analysis software.

MassHunter acquisition software has two possible configurations:

- Workstation (standalone system with no content management) or,
- Networked Workstation (client computer with content management on a separate server)
- Prepare the MassHunter Workstation computer the same way, regardless of the configuration.

Installation

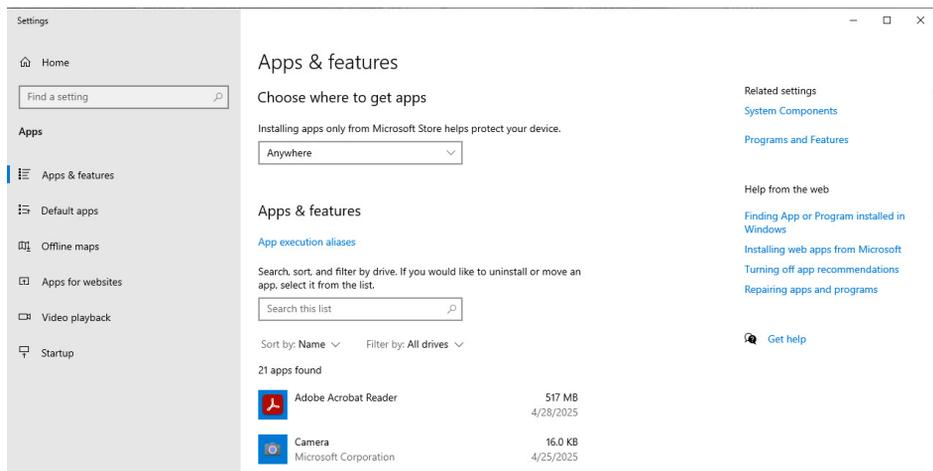
Step 1: Prepare for Installation

Step 1: Prepare for Installation

- 1 Run **Windows Update** to ensure that the latest critical updates and security fixes are installed. **Windows Update** must be complete before installing MassHunter Workstation Qualitative or Quantitative Analysis.



- 2 If upgrading MassHunter Workstation Quantitative Analysis B.05.00 or earlier, use the Uninstall Programs command in the Windows Control Panel to remove the programs to be upgraded. See Microsoft Windows help for more information on uninstalling programs.



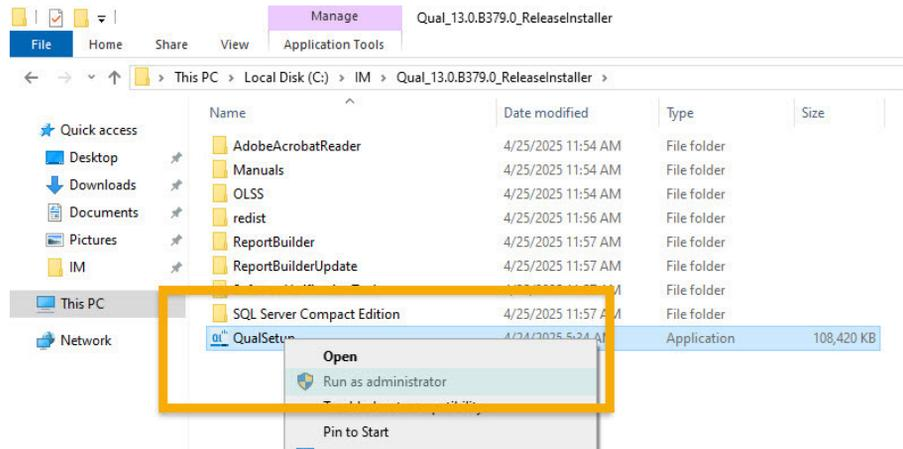
- 3 Once the programs are removed, reboot the computer.

Installation

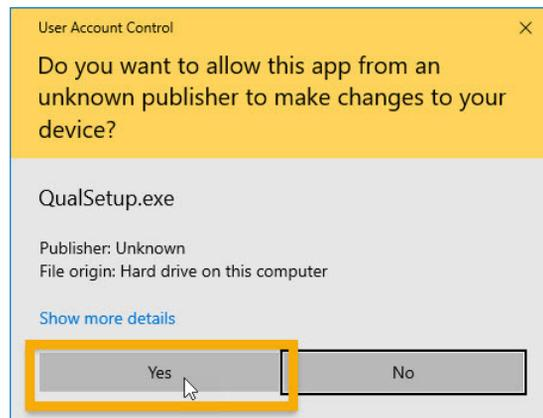
Step 2: Install the Qualitative Analysis Program

Step 2: Install the Qualitative Analysis Program

- 1 From the installation media, right-click QualSetup.exe and select **Run as administrator**.



- 2 Click **Yes** to *Do you want to allow this app to from an unknown publisher to make changes to your device?*

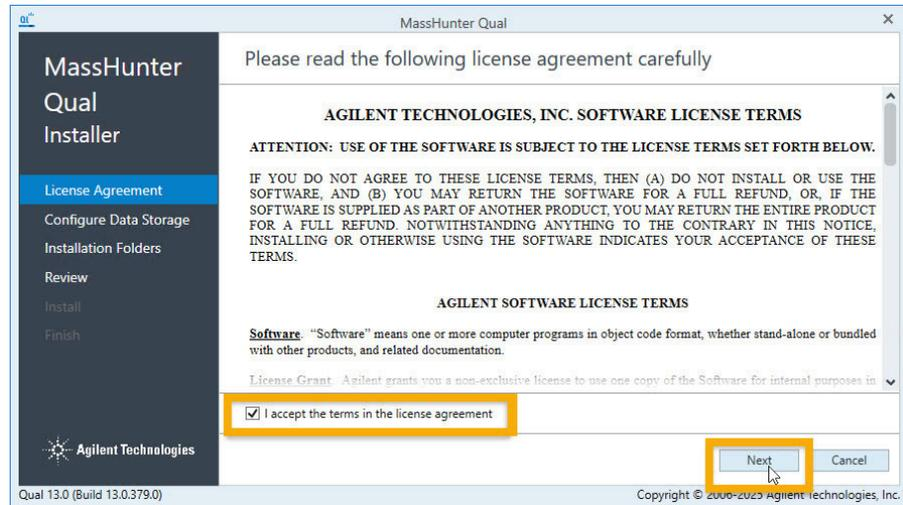


Installation

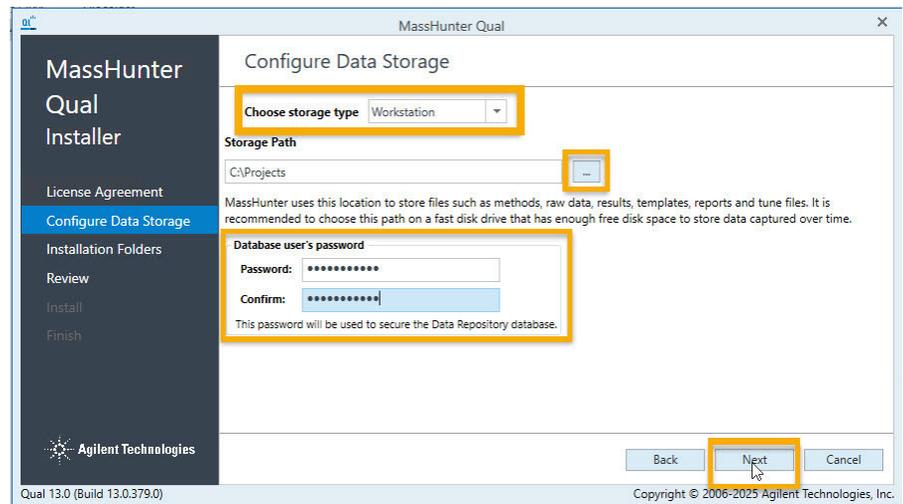
Step 2: Install the Qualitative Analysis Program

- 3 Click the check box to accept the terms of the license agreement, then click **Next**.

Through the installation, click **Next** to continue with the installation, **Back** to review previous selections, or **Cancel** to end the installation.



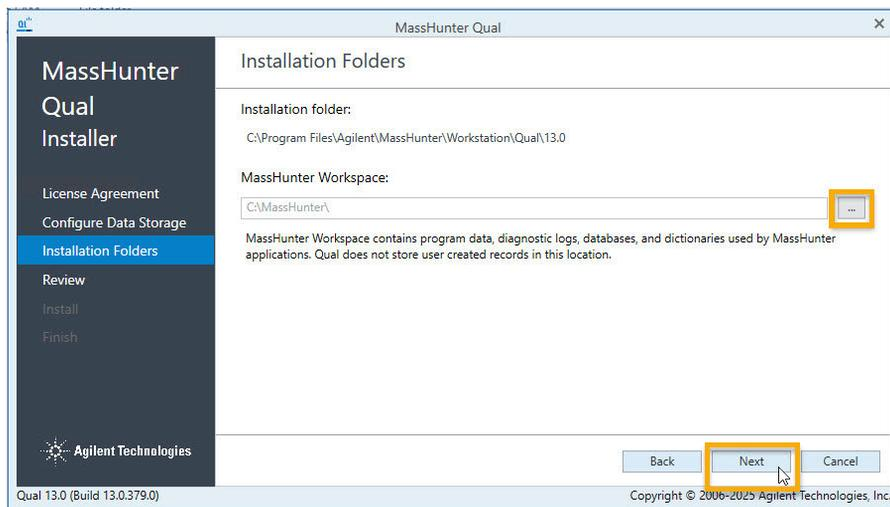
- 4 Configure the desired storage type, path and enter a password. Confirm the password and click **Next**.



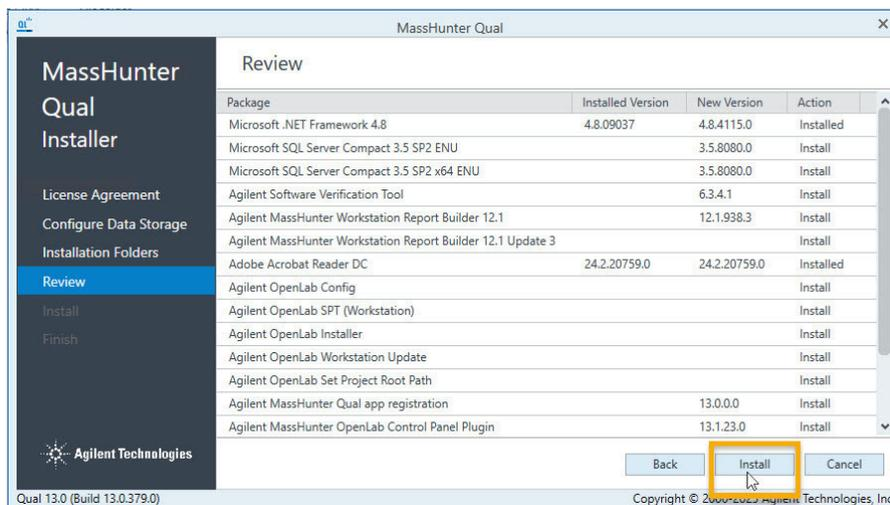
Installation

Step 2: Install the Qualitative Analysis Program

- 5 Use the default location for the software files or click  to choose a different location, then click **Next**.



- 6 Review the installation information. Click **Back** to change any settings, or click **Install** to continue.
Installation may take a few minutes.

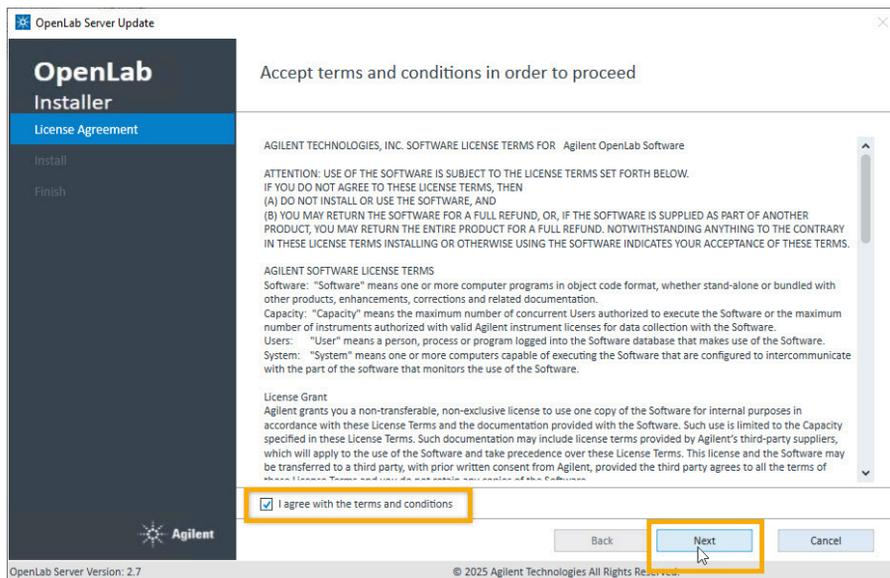


Installation

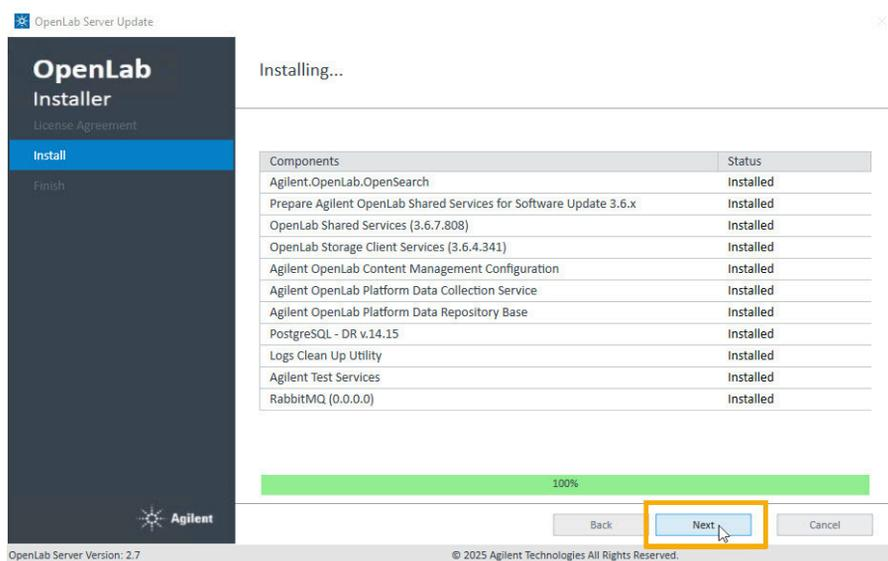
Step 2: Install the Qualitative Analysis Program

- 7 Click the check box to accept the terms of the OpenLab license agreement, then click **Next**.

Installation may take a few minutes.



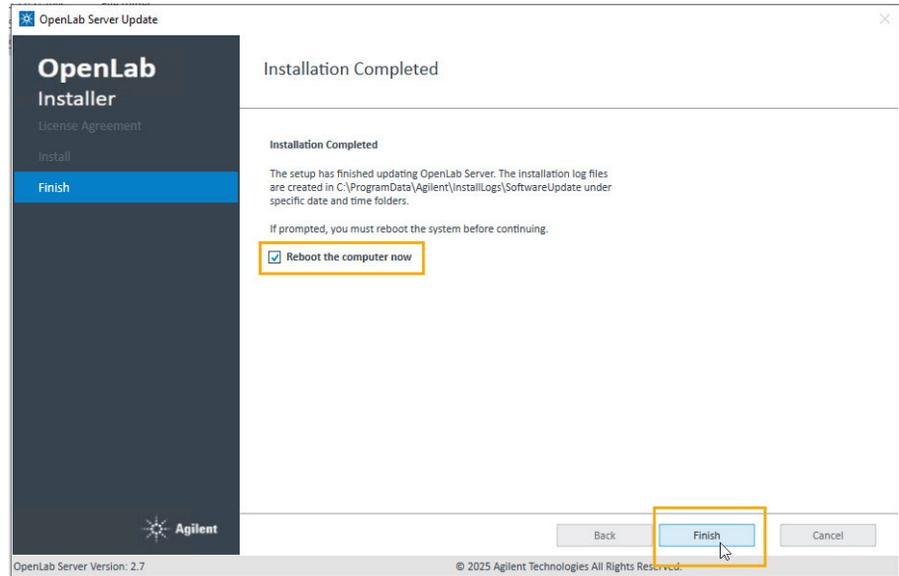
- 8 Click **Next** when installation is complete.



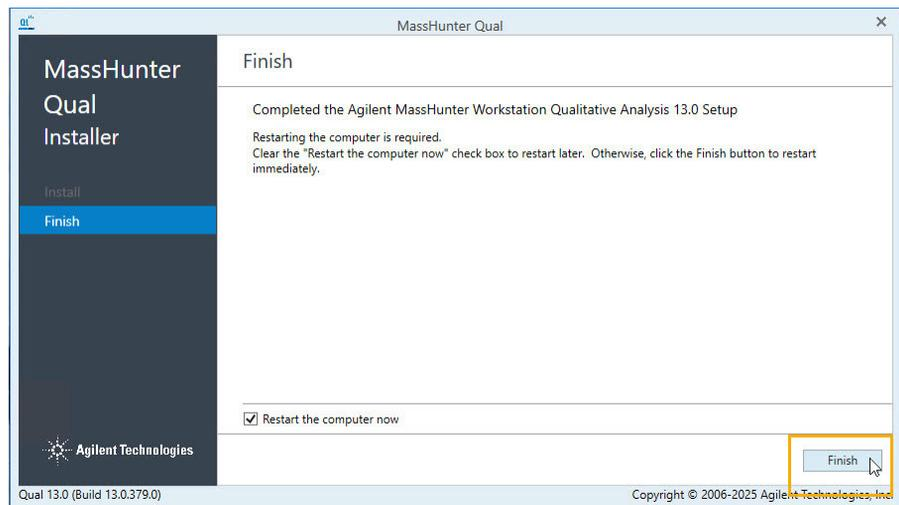
Installation

Step 2: Install the Qualitative Analysis Program

- 9 Click **Finish**. The computer will reboot.



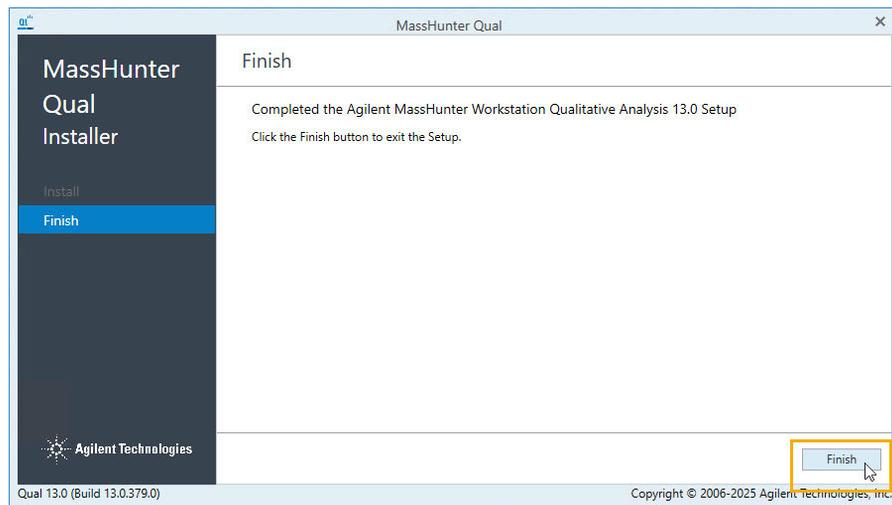
- 10 Upon login, the installation will continue automatically. When completed, click **Finish**. The computer will reboot.



Installation

Step 2: Install the Qualitative Analysis Program

- 11 Upon login, the installation will continue automatically. When completed, click **Finish**.

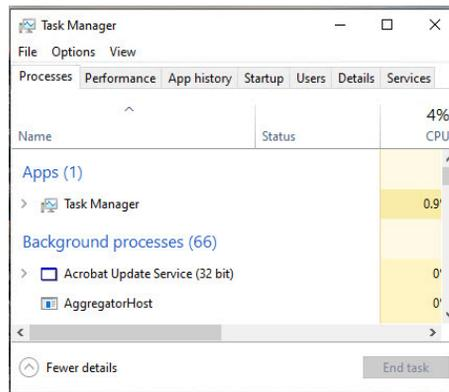


Installation

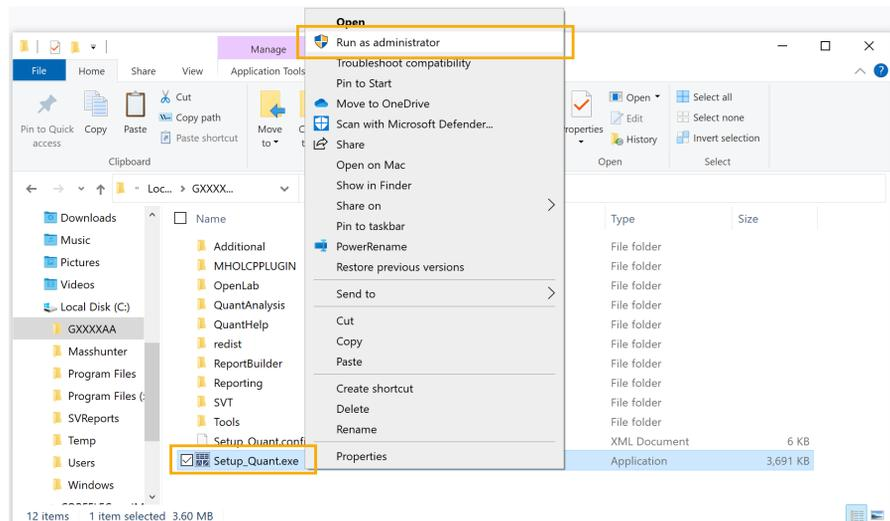
Step 3: Install or Upgrade the Quantitative Analysis Program

Step 3: Install or Upgrade the Quantitative Analysis Program

- 1 Verify that no other programs are running the system, including Windows Update.



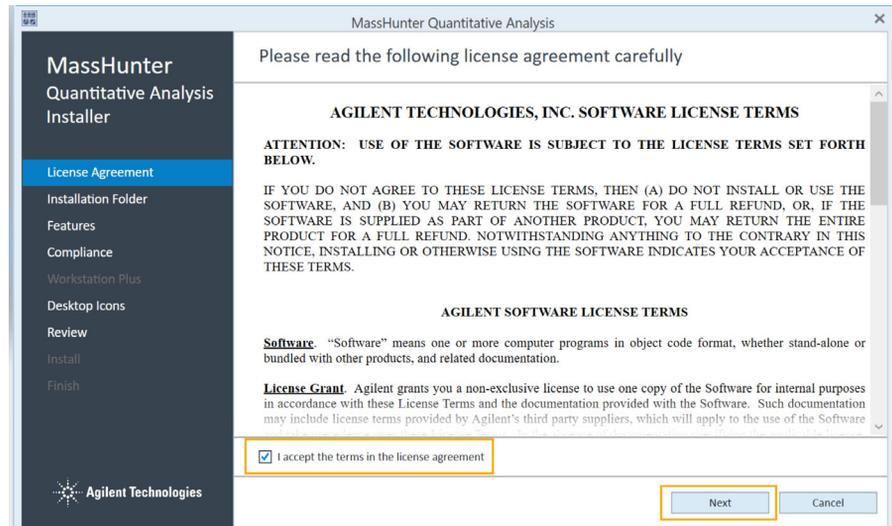
- 2 From the installation media, right-click **Setup_Quant.exe**, and select **Run as administrator**.



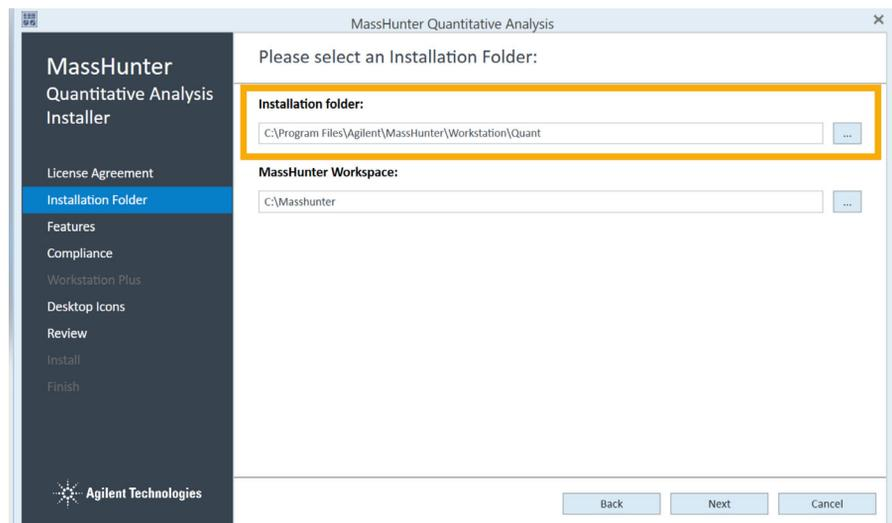
Installation

Step 3: Install or Upgrade the Quantitative Analysis Program

- 3 Review the License Agreement, and accept the terms. Click **Next** to continue.



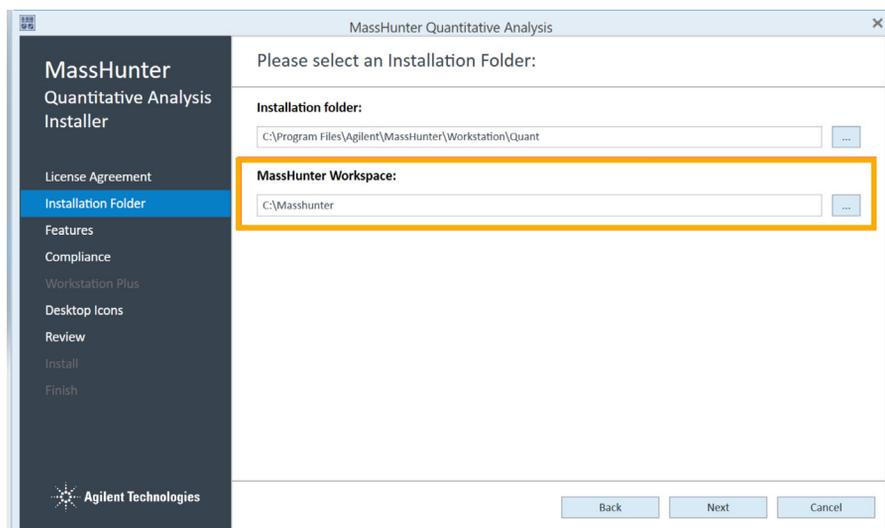
- 4 Install the program to the default location, or click **Browse** to choose a different location.



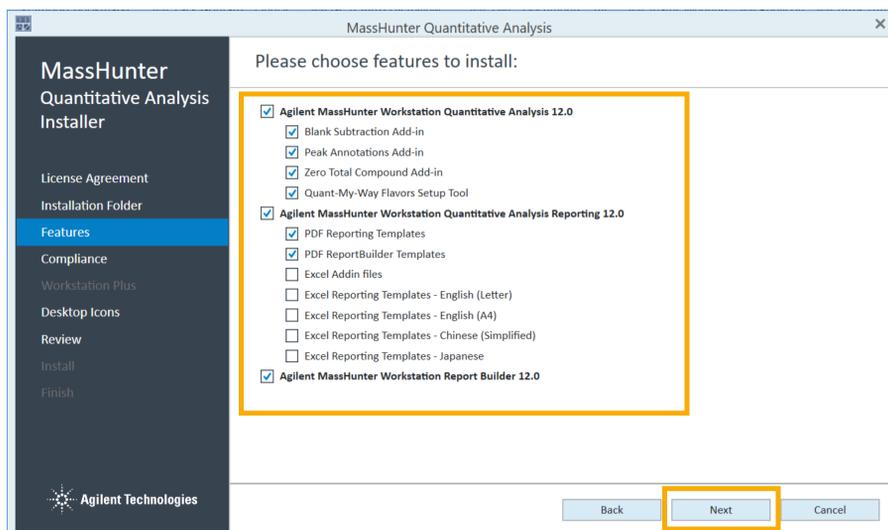
Installation

Step 3: Install or Upgrade the Quantitative Analysis Program

- 5 The MassHunter workspace installs to the default location, a shared directory utilized by multiple MassHunter products. Changing this location is not recommended if other MassHunter applications are already installed. Click **Next** to continue.



- 6 Select the features to install, and click **Next** to continue.



Installation

Step 3: Install or Upgrade the Quantitative Analysis Program

Use Add-ins for additional features:

- **Blank Subtraction Add-in:** This add-in allows subtraction of the 'Blank' and/or 'DoubleBlank' sample final concentration and places the blank-subtracted concentrations into the target compound Custom Calculation column of the batch table. If a batch contains multiple 'Blank' and/or 'DoubleBlank' samples, the average blank concentration is used.
- **Peak Annotations Add-in:** In the Quantitative Analysis program, this add-in allows an abbreviation to be added to the batch table that indicates the reason for manually integrating target or ISTD compounds.
- **Zero Total Compound Add-in:** In the Quantitative Analysis program, this add-in allows the zeroing out of the peaks for all compounds in a group.
- **Quant-My-Way Flavors Setup Tool:** This tool allows use of the Quant-My-Way Flavors Setup program to create additional Quantitative Analysis flavors that display the tools and menu options needed for the laboratory or function. See the Agilent MassHunter Quantitative Analysis Quant-My-Way Flavors Setup Familiarization Guide.

The reporting templates are installed on the local hard drive with the program to MassHunter/Report Templates/Quant/. If planning to use MassHunter Compliance Software, these templates must be uploaded to ECM before they can be accessed from the Quantitative Analysis program.

Installation

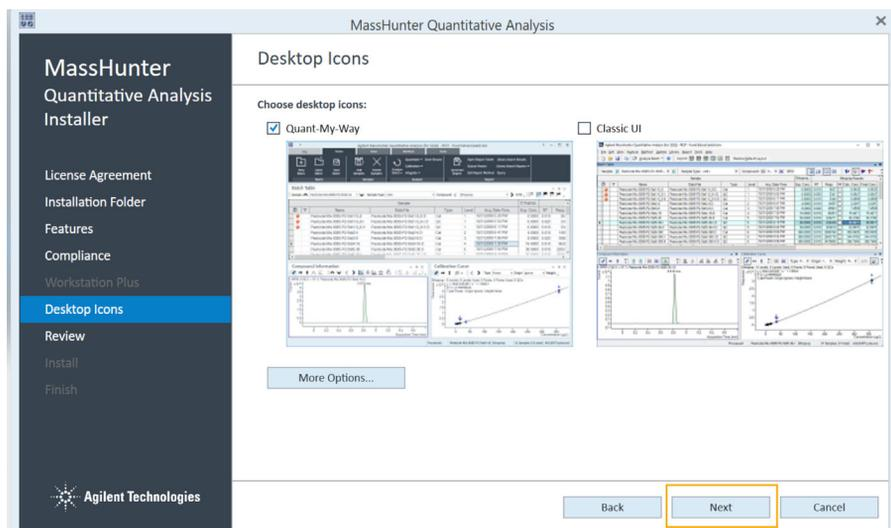
Step 3: Install or Upgrade the Quantitative Analysis Program

- 7 Select the desktop icons, and click **Next** to continue.

The default set of icons open Quantitative Analysis in the **Quant-My-Way** user interface and include these standard flavors:

- Quantitative Analysis MS
- Quantitative Analysis QQQ
- Quantitative Analysis TOF
- Quantitative Analysis Q-TOF
- Unknowns Analysis
- Vanilla Pre-Configured Flavor

To install the standard set of icons for the **Classic** interface, select **Classic UI**. Choose to install a single set of interface icons, or select both interfaces to install.

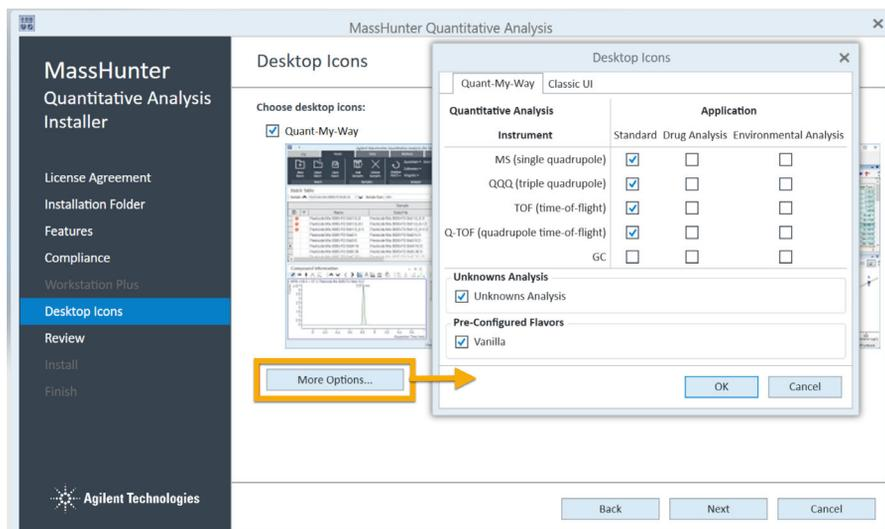


To include a GC instrument user interface or to manually select which desktop icons to install, click **More Options**, and select each desktop icon to install.

Installation

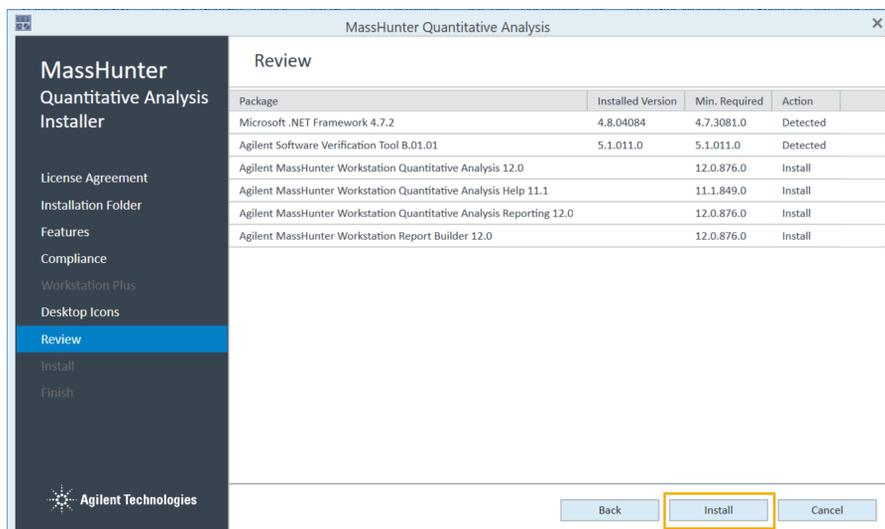
Step 3: Install or Upgrade the Quantitative Analysis Program

The **Quant-My-Way** user interface offers a pre-configured **Vanilla** flavor. This flavor is based on the standard QQQ flavor and provides a simplified user interface that includes only the features needed for basic data analysis.



After installation, change which icons are displayed. See the Online Help for more information.

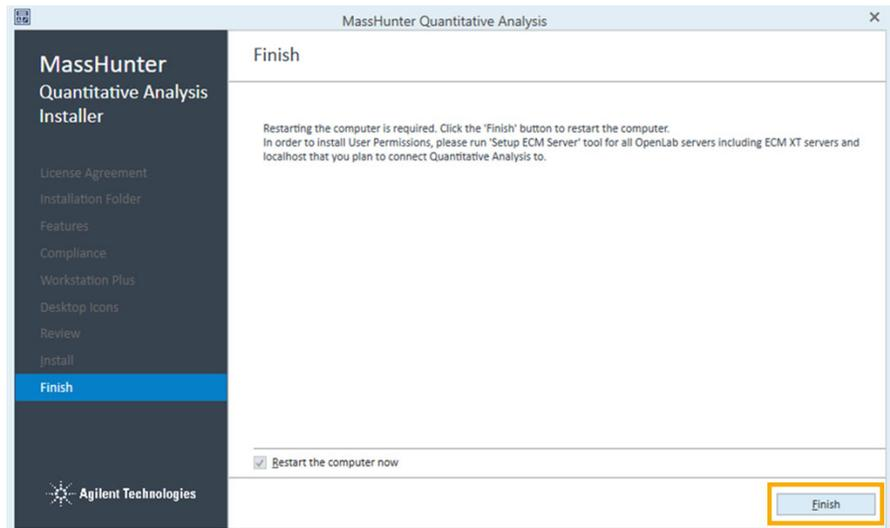
- 8 Click **Back** to change the settings, or click **Install** to continue. Installation may take a few minutes.



Installation

Step 3: Install or Upgrade the Quantitative Analysis Program

9 Click **Finish**.



10 If the **Service Packs** folder exists on the installation media:

- a Open the **Service Packs** folder. If the folder contains another folder, such as SP1, open that folder.
- b Double-click the installation file ending with **SPx.msi**.
- c Follow the instructions to install the service pack.

11 To complete the installation, restart the machine.

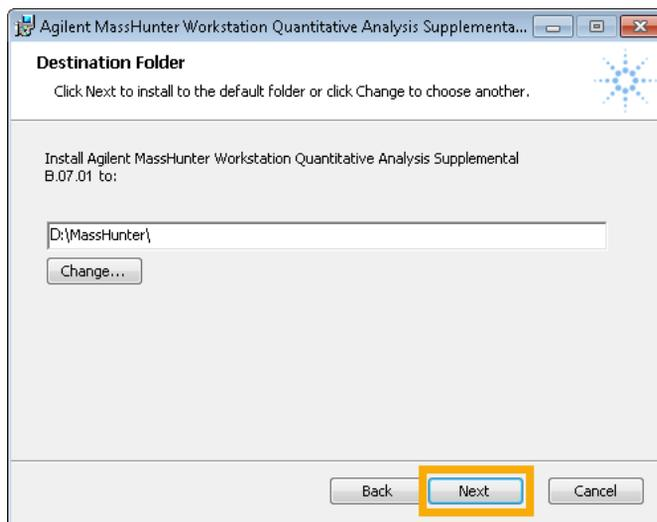
Installation

Step 4: Install Quantitative Analysis Supplemental Items

Step 4: Install Quantitative Analysis Supplemental Items

Perform this step to copy the supplemental manuals, report template index, technical videos, and data to the computer from the installation media.

- 1 From the Quantitative Analysis installation media, right-click **Supplemental.msi**, and select **Install**.
- 2 Review the terms, and select **I accept the terms in the License Agreement**.
- 3 Click **Advanced** or **Install**
 - Click **Install** to continue the installation using the default options for all users.
 - Click **Advanced** to change the installation options.
 - a On the **Destination Folder** page, the default location is **D:\MassHunter**. Click **Change** to choose a different location for the supplemental items. Click **Next** when the location is set.

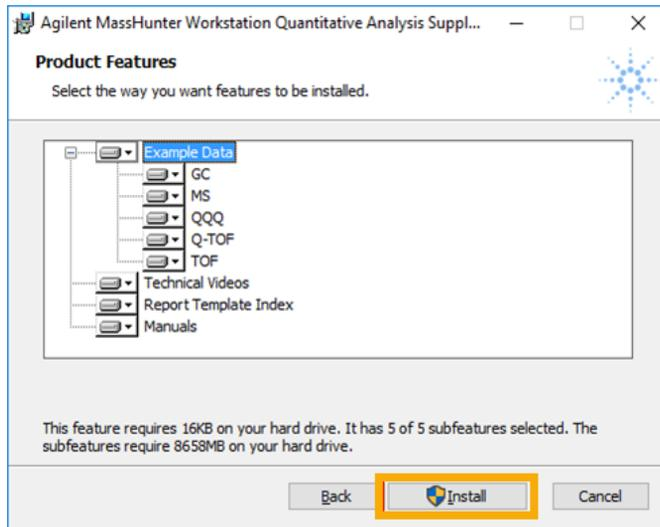


- b On the **Product Features** page, select the way to install each feature.

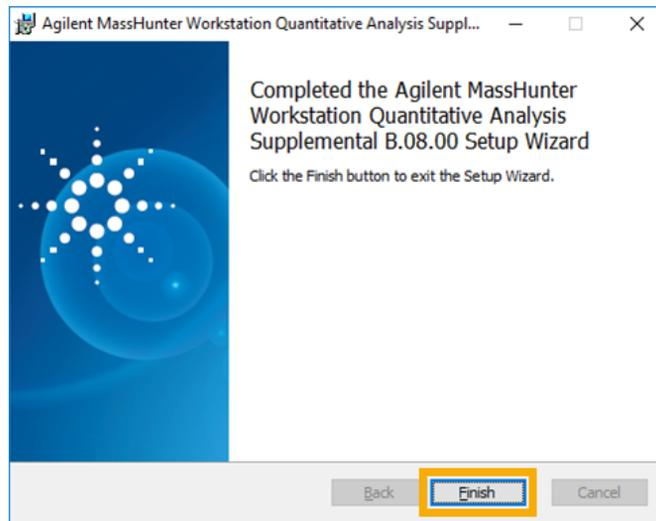
Installation

Step 4: Install Quantitative Analysis Supplemental Items

c Click **Install**.



4 Click **Finish**.



To view the supplemental materials:

- For Windows 7, select **Start > All Programs > Agilent > MassHunter Workstation > Quantitative Analysis Supplemental**.

Installation

Step 4: Install Quantitative Analysis Supplemental Items

- For Windows 10 and 11, select **Start > All apps > Agilent MassHunter Quantitative > Quantitative Analysis Supplemental**.

Installation

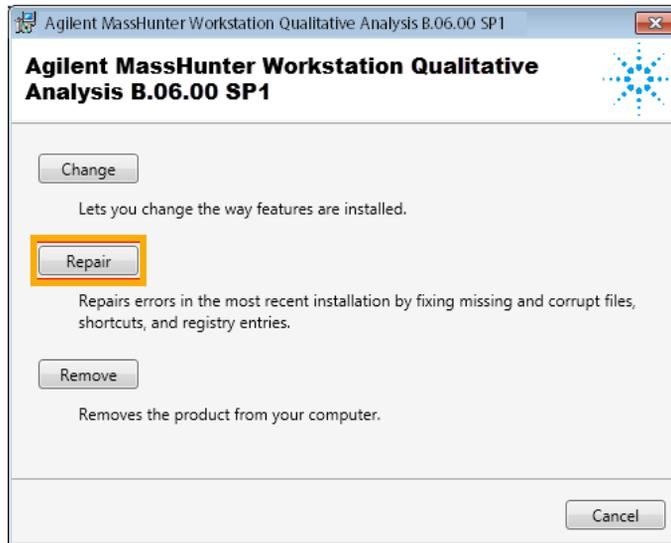
Step 5: Configure Excel for MassHunter

Step 5: Configure Excel for MassHunter

Excel for Qualitative Analysis

Do this step only if Microsoft Excel was upgraded after Qualitative Analysis or Quantitative Analysis was installed or upgraded.

- 1 Click **Repair**.



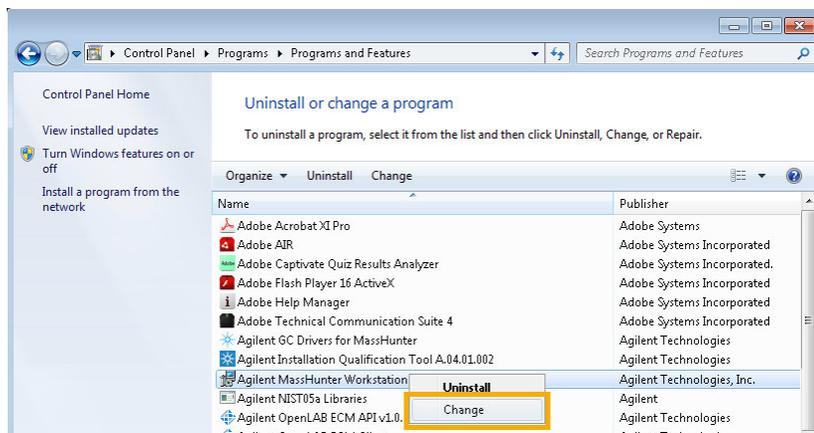
- 2 Follow the instructions to complete the process.

Configure Excel for Quantitative Analysis

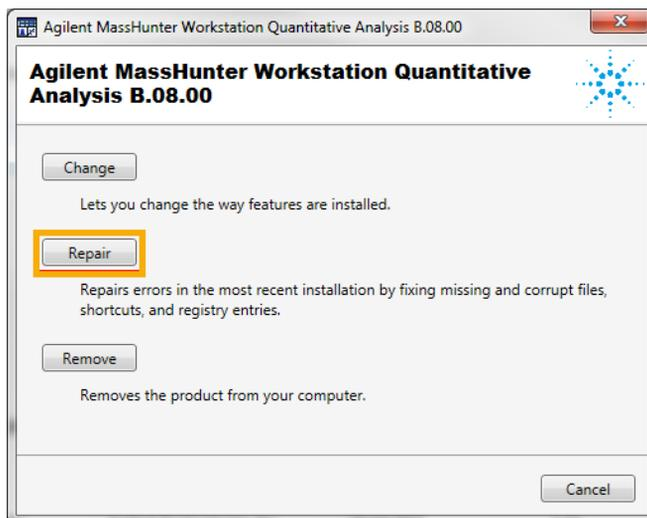
- 1 Click **Start > Control Panel > Programs and Features**.
- 2 Right-click **Agilent MassHunter Workstation Quantitative Analysis**, and select **Change**.

Installation

Step 5: Configure Excel for MassHunter



3 Click **Repair**.



4 Follow the instructions to complete the process.

www.agilent.com

© Agilent Technologies, Inc. 2025

D0130987

May 2025

