



MassHunter Explorer

Installation Guide

Notices

Document Identification

D0109800
February 2024 A.01

Copyright

© Agilent Technologies, Inc. 2024

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This guide is valid for the 1.0 revision or higher of MassHunter Explorer until superseded.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.

History

Revision	Date	Change
A.01	February 2024	Initial release.
A.01	February 2024	Corrected hyperlinks.

In This Guide

This guide provides licensing and installation instructions for MassHunter Explorer.

Contents

- 1 Licensing and Download 7**
 - About MassHunter Explorer Software Licensing 8**
 - Software Maintenance Agreement (SMA) 8
 - Requirements 9**
 - PC Requirements 10**
 - Minimum PC Requirements 10
 - Setting up an Agilent.com account 11
 - Download the software 13
- 2 Installation 17**
 - Install MassHunter Explorer Software 18
 - Activate MassHunter Explorer Software 27
 - Offline Activation 29
- 3 Getting Started with MassHunter Explorer Software 33**
 - Next Steps 34**
 - User Documentation 34
 - Agilent Community 34
- 4 Transferring MassHunter Licenses 35**
 - Returning MassHunter Explorer software licenses and uninstalling 35

This page intentionally left blank.



1

Licensing and Download

This chapter describes how to obtain and activate a MassHunter Explorer software license and download the installation files.

About MassHunter Explorer Software Licensing

Software Maintenance Agreement (SMA)

Bundled into MassHunter Explorer software is a one-year software maintenance agreement.

The software maintenance agreement does not include installation services. Installation or upgrade services must be purchased by contacting your sales representative in your region.

Requirements

To receive a license, you must:

- Be named as the software, install, or contact person during purchase.
- Have an Agilent.com account.
- Obtain the entitlement certificate ID for this software from the License and Delivery portal.

Before you begin:

- Ports 7070 and 7071 need to be open for licensing to work. If other software use the same ports, it needs to be uninstalled before activation of the MassHunter Explorer software.
- If you **do not** have an Agilent.com account, an email invitation to set a password is sent to the individual indicated as the point of contact after purchase. The email address is no-reply@mailing.agilent.com with the title *Agilent Software and Licensing Portal: Account and Password Verification*.
- If you **do have** an account, an email with the Entitlement Certificate is automatically sent to the individual indicated as the point of contact. This email address is donotreply_agilent@flexnetoperations.com with the title *Your Agilent Entitlement Information*.

If you have not received the expected email for your product, contact your vendor or internal support.

PC Requirements

Minimum PC Requirements

The following are the minimum computer hardware, software, and network requirements to run Agilent MassHunter Explorer software.

Item	
Description	High Capacity MassHunter-ready Computer (M6082AA)/Hewlett-Packard Z4 G4 Minitower
Processor speed (CPU)	Intel Xeon W-2235, 6 core, 3.8 GHz
Physical memory (RAM)*	64 GB
Hard Disk	1 TB M.2 NVMe SSD - Primary (C:\) Boot. 6 TB x 4 RAID10 (12 TB) - Data (D:\)
Graphic Resolution	1920 x 1080
Graphics Card	NVIDIA Quadro T400 4 GB (3)mDP GFX w/
Operating System	Windows 10 or 11 Pro

* Allocate 4GB of RAM per logical thread to ensure that the system has sufficient total RAM to accommodate the computational workload.

Setting up an Agilent.com account

An Internet connection is needed to access Agilent.com.

NOTE

If you are an existing customer with an Agilent.com account, skip to “**Download the software**” on page 13.

- 1 After purchase, if you are a new customer, an email is sent with instructions to verify and set a password for an Agilent.com account from **no-reply@mailing.agilent.com**. Click the link to verify the email account and set a password.



Welcome to the Agilent Software and Licensing Portal! Your account has been created either through a recent purchase or request to be added to our portal.

Please click on [this link](#) to verify your account and set a password.

If you have software or licensing available to download and activate, you will receive a separate email detailing this information.

For questions or comments, please contact us:

[Contact Us | Agilent](#)

Thank You,
Your Agilent Team

NOTE

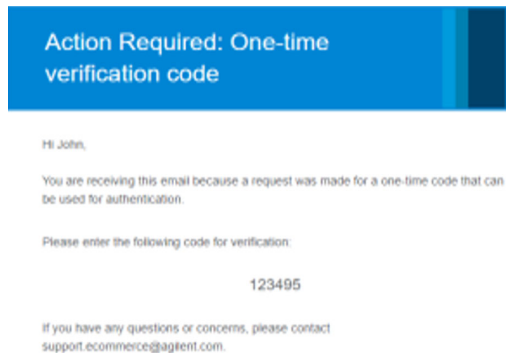
If you do not receive an email, check your email spam and junk folders. If the email is not found, contact the Agilent Global Service Contact Center (<https://www.agilent.com/en/contact-us/page>).

Licensing and Download

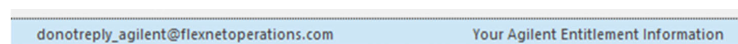
Setting up an Agilent.com account

- The Create a Password page loads in a browser window. Enter a password and confirm. Click **Create Account**.

- Once the password is set, an email is sent with a one-time verification code to verify the account. Enter the code to complete the verification to Agilent.com.



An email is sent with the Entitlement Certificate to the verified email address from **donotreply_agilent@flexnetoperations.com** with the subject line *Your Agilent Entitlement Information*.



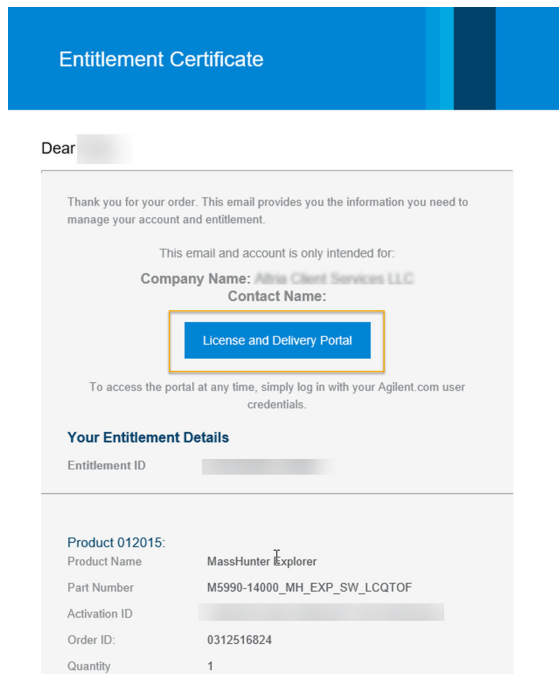
NOTE

If you do not receive an email, check your email Spam and Junk folders. If the email is not found, please reach out to your local Agilent sales and support contact.

Download the software

An Internet connection is needed to access the License and Delivery Portal.

- 1 From within the Entitlement email, click the **License and Delivery Portal** button.

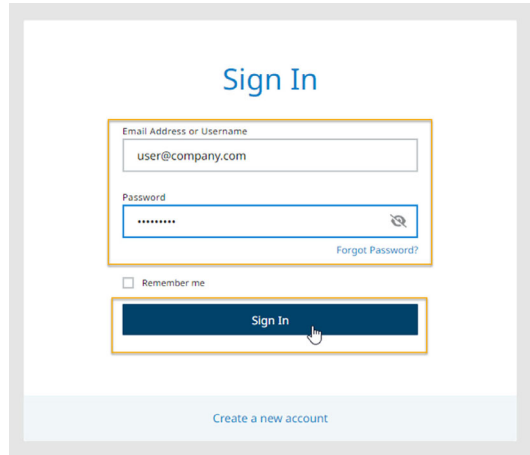


1

Licensing and Download

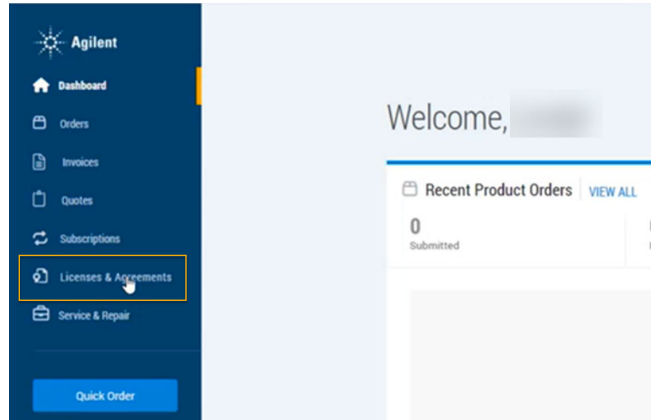
Download the software

- Using the Sign In screen, enter your login credentials and click **Sign In** to log into Agilent.com to access the License and Delivery portal.



The image shows a 'Sign In' web form. At the top, the text 'Sign In' is displayed in blue. Below it, there are two input fields: 'Email Address or Username' containing 'user@company.com' and 'Password' containing several asterisks. A 'Forgot Password?' link is located to the right of the password field. Below the password field is a checkbox labeled 'Remember me'. At the bottom of the form is a dark blue button with the text 'Sign In' and a mouse cursor icon. Below the main form area, there is a link that says 'Create a new account'.

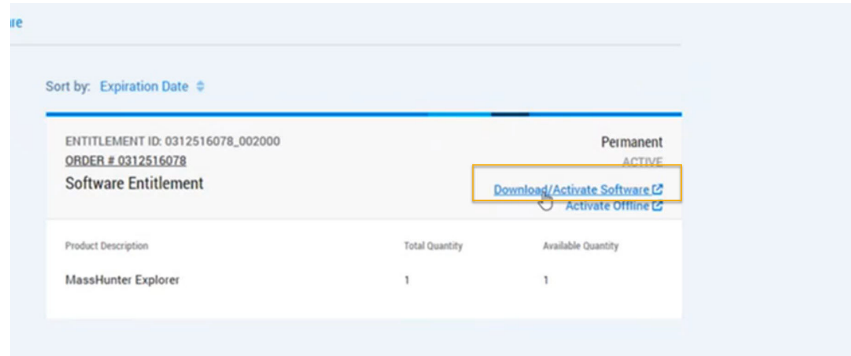
- Once in the portal, click **Licenses & Agreements**.



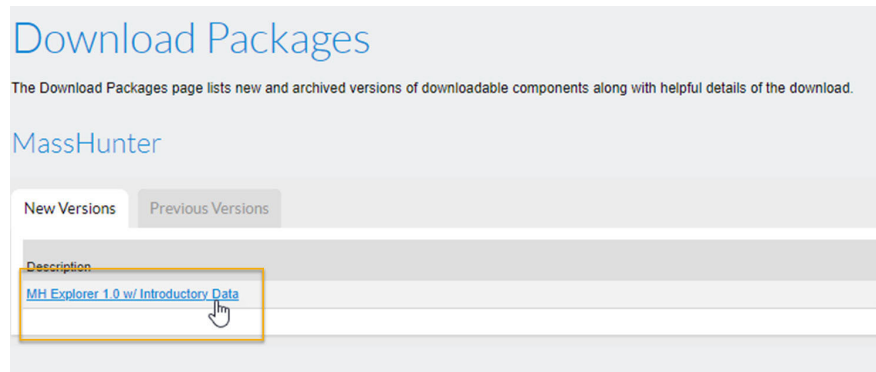
Licensing and Download

Download the software

- 4 Under Software Licenses, locate the MassHunter Explorer software listing, then click **Download/Activate Software**.



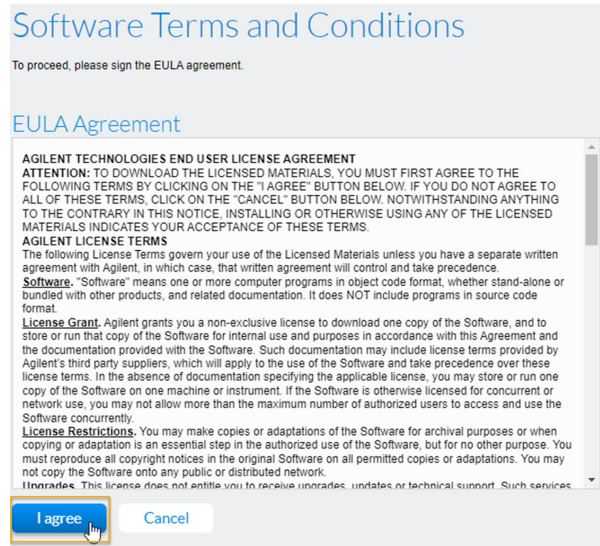
- 5 The Download Packages screen loads. Click the **MH Explorer 1.0 w/Introductory Data** link.



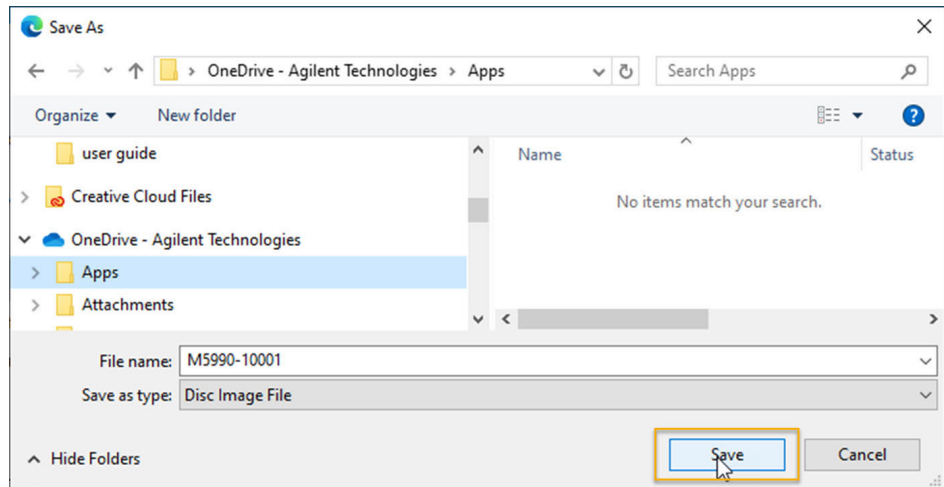
Licensing and Download

Download the software

- 6 Accept the Software Terms and Conditions by clicking **I Agree**.



- 7 In the downloads notification of the browser, click **Save As**, browse to a known location on the installation computer, then click **Save**.



You are now ready to install the software.



2 Installation

This chapter describes how to install the MassHunter Explorer software.

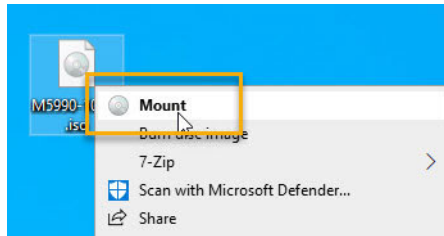
Install MassHunter Explorer Software

WARNING

These instructions should only be used by customers with standalone workstations. MassHunter Explorer cannot be installed on networked workstations.

Before installing the MassHunter Explorer software, you must uninstall any previous version. See [“Returning MassHunter Explorer software licenses and uninstalling”](#) on page 35.

- 1 Navigate to the directory where the MassHunter Explorer software installation file is saved. Right-click, then select **Mount**.



This action opens a virtual drive that is visible in File Explorer.

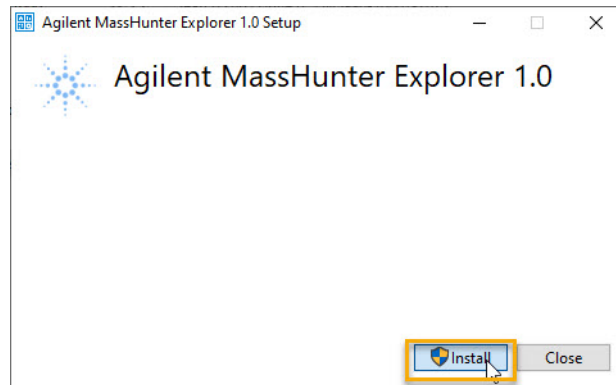
- 2 Select all the files in the mounted drive, then copy and paste them to a directory on the C:\ drive.
- 3 From the directory on the C:\ drive, double click to run the setup program, **MassHunterExplorerSetup.exe**.

Name	Date modified	Type	Size
Support	15/12/2023 6:53 AM	File folder	
Data - Explorer Basic Introduction.zip	14/12/2023 8:26 PM	Compressed (zipp...	1,882,958 KB
LCMS User Documentation Information...	1/12/2023 11:28 PM	Microsoft Edge P...	522 KB
MassHunterExplorerSetup.exe	12/12/2023 7:52 AM	Application	698,592 KB

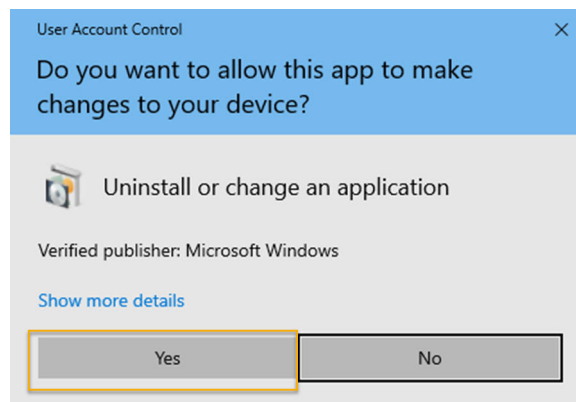
Installation

Install MassHunter Explorer Software

- 4 Click **Install**.



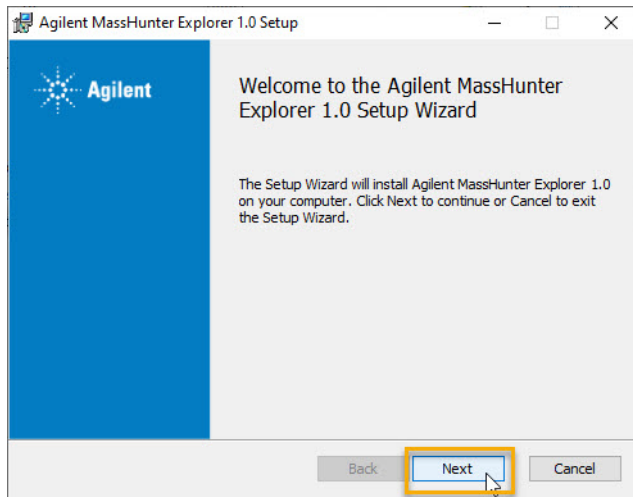
- 5 Click **Yes** to *Do you want to allow this app to make changes to your device?*



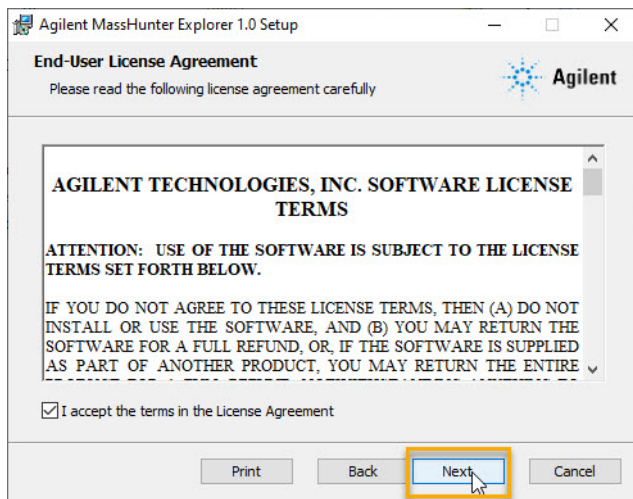
Installation

Install MassHunter Explorer Software

- 6 Click **Next**.



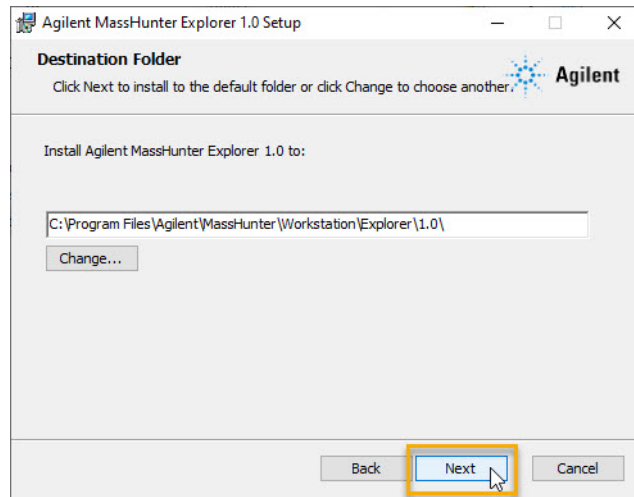
- 7 Click the check box to accept the terms of the license agreement, then click **Next**.



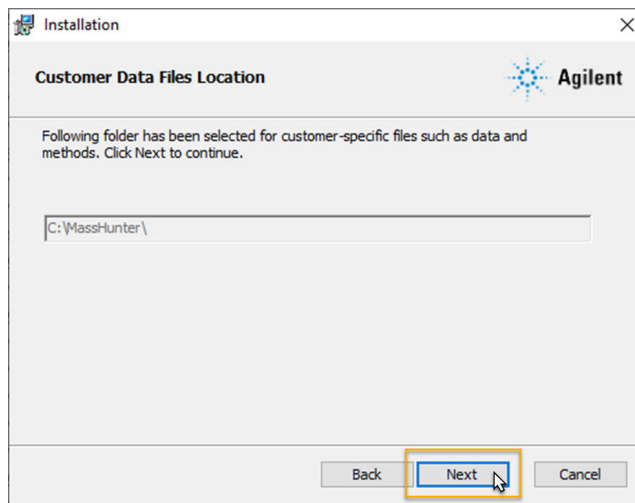
Installation

Install MassHunter Explorer Software

8 Click **Next**.



9 Click **Next**.

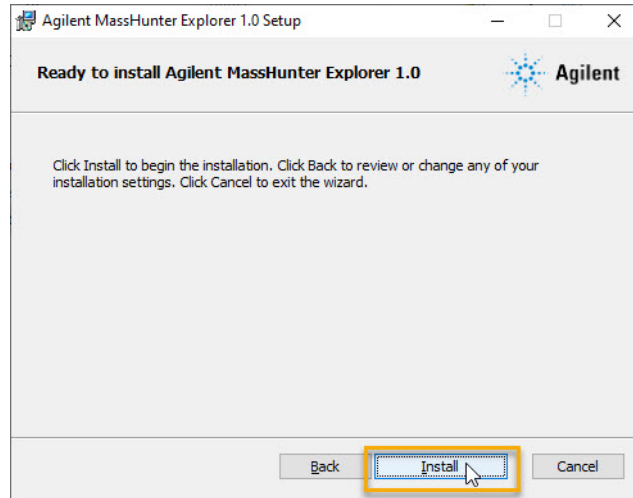


Installation

Install MassHunter Explorer Software

10 Click **Install**.

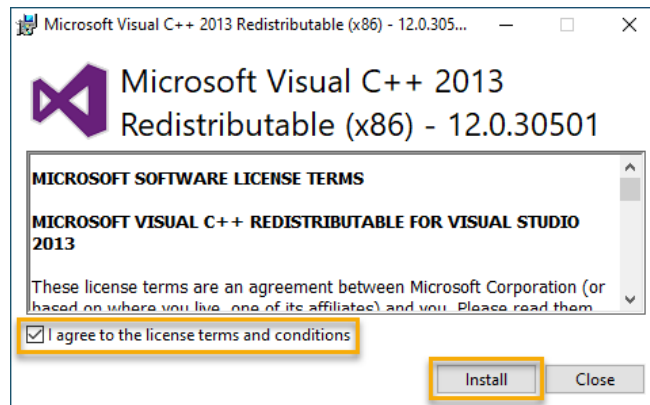
When you click **Install**, the installation of the required packages will start.



NOTE

If .NET framework or Agilent OpenLab licensing is not installed on your system, its installation is automatically triggered by the installation wizard. Skip to step 18 if already installed.

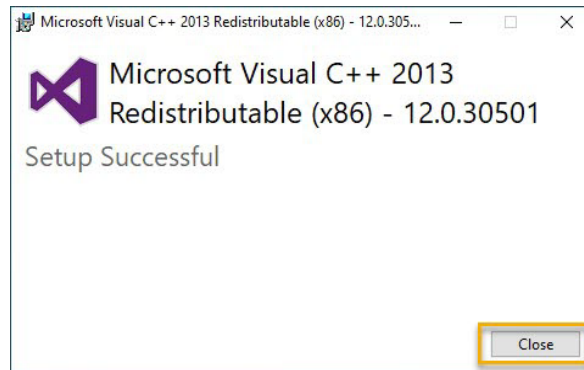
11 Click to agree to the license terms and conditions, then click **Install**.



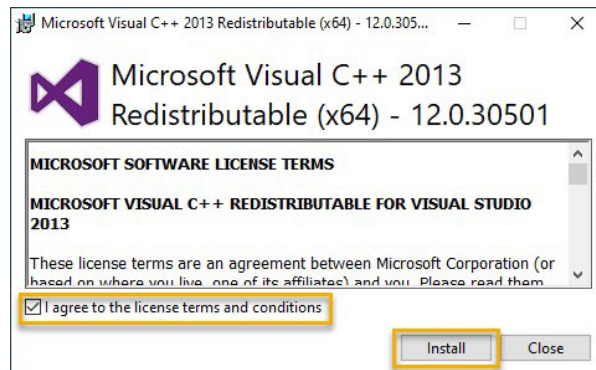
Installation

Install MassHunter Explorer Software

12 Click **Close**.



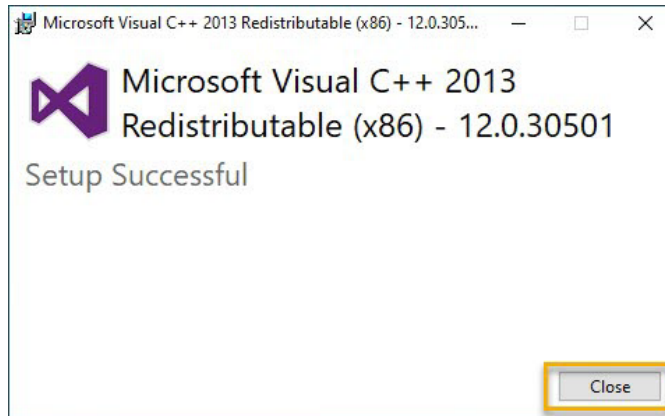
13 Click to agree to the license terms and conditions, then click **Install**.



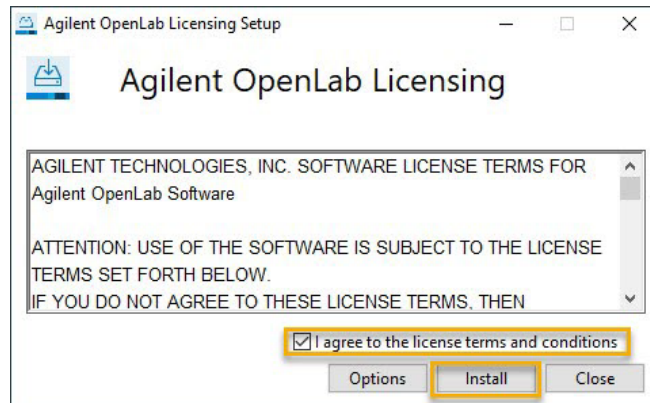
Installation

Install MassHunter Explorer Software

14 Click **Close**.



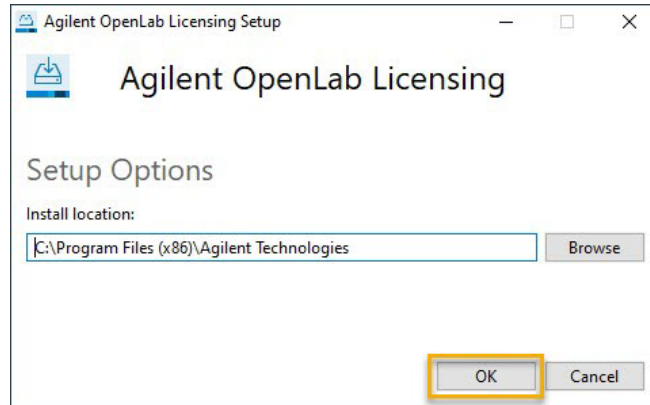
15 Click to agree to the license terms and conditions, then click **Install**.



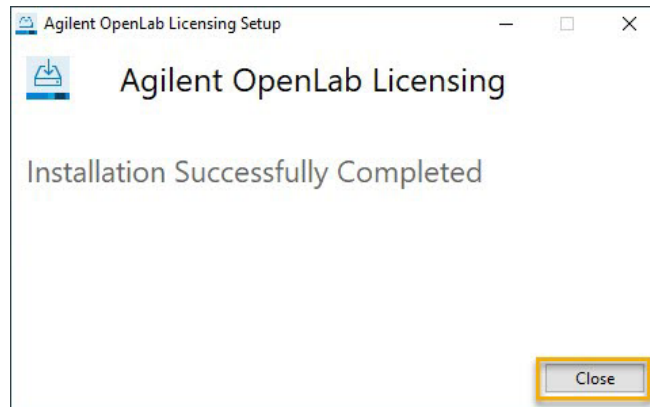
Installation

Install MassHunter Explorer Software

16 Click **OK**.



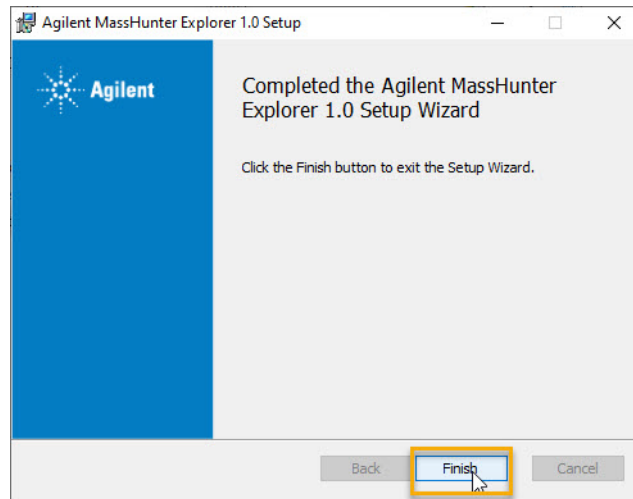
17 Click **Close**.



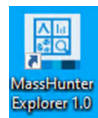
Installation

Install MassHunter Explorer Software

18 Click **Finish**.



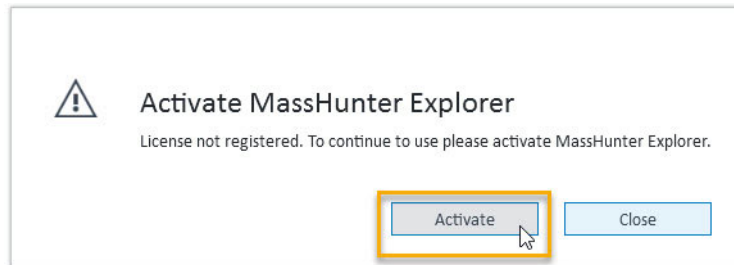
19 When the installation is complete, the MassHunter Explorer icon is displayed on your desktop.



Activate MassHunter Explorer Software

The installation PC must be connected to a network with Internet access to activate the license. If the installation PC is not Internet connected, see **“Offline Activation”** on page 29.

- 1 Double click the **MassHunter Explorer 1.0** icon to open the program.
- 2 Upon launch, the Activate MassHunter Explorer dialog box displays. Click **Activate**. The Agilent Licensing window displays in a browser window.



- 3 Verify that the type of License Activation method is **Online Activation**.
- 4 From the entitlement email, copy the **Activation ID**.

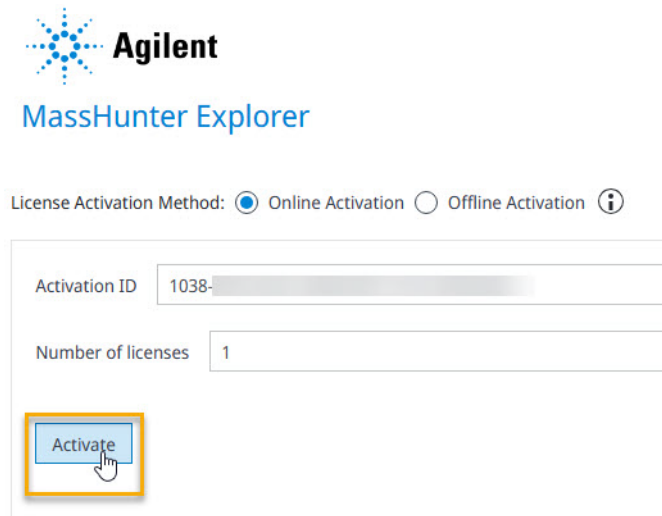
Product 012015:	
Product Name	MassHunter Explorer
Part Number	M5990-14000_MH_EXP_SW_LCQTOF
Activation ID	1038- eec
Order ID:	0312516824
Quantity	1
Permanent	Yes

- 5 Paste the Activation ID into the **Activation ID** field.
- 6 Ensure that the License Activation option is selected.

Installation

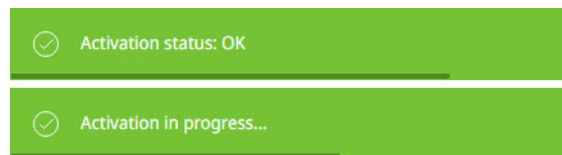
Activate MassHunter Explorer Software

- 7 Enter the number of licenses to be used, and click **Activate**.



The screenshot shows the Agilent MassHunter Explorer activation window. At the top is the Agilent logo and the text "MassHunter Explorer". Below this, the "License Activation Method" is set to "Online Activation" (indicated by a selected radio button) with an information icon to its right. The "Offline Activation" option is unselected. There are two input fields: "Activation ID" containing "1038-" and "Number of licenses" containing "1". A blue "Activate" button is highlighted with a yellow border and a mouse cursor pointing to it.

A confirmation message will flash along the bottom of the screen.



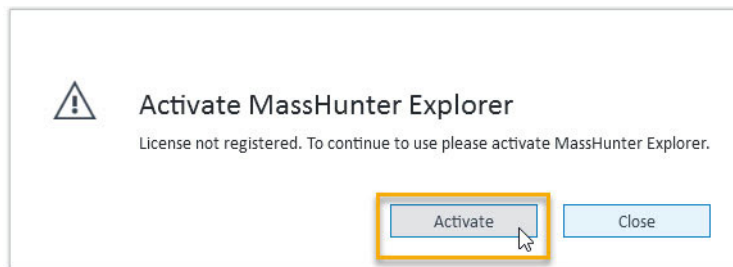
- 8 Once complete, close the browser window.

The MassHunter Explorer software is now licensed and ready to use.

Offline Activation

Use this activation procedure if your workstation is not connected to the Internet.

- 1 Double click the **MassHunter Explorer 1.0** icon to open the program. Upon launch, the Activate MassHunter Explorer dialog box displays.
- 2 Click **Activate**. The Agilent Licensing window displays in a browser window.



- 3 Select the License Activation method **Offline Activation**.

License Activation Method: Online Activation Offline Activation ⓘ

- 4 From the entitlement email, copy the **Activation ID**.

Product 012015:	
Product Name	MassHunter Explorer
Part Number	M5990-14000_MH_EXP_SW_LCQTOF
Activation ID	1038- eec
Order ID:	0312516824
Quantity	1
Permanent	Yes

5 Add the Activation ID into the **Activation ID** field.

Generate Activation Request | Install Activation Response | Activation Confirmation

Step 1 Generate Activation Request file:

- Enter Activation ID's and quantities for all licenses you want to activate offline. Click the + button to add these to the table.
- Update existing Activation ID quantities in the table by editing the value in the Total column.
- When all Activation ID's and quantities are finalized, click **Generate Activation Request** button to create an *Activation Request* file.

Activation ID	Number of licenses
ad1b-	1

Generate Activation Request

6 Enter the **Number of licenses**.

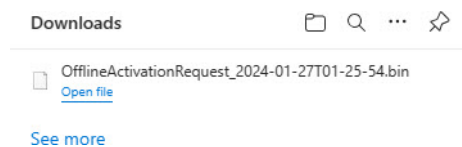
7 Click the + button to display the entered information in the license table. Update existing Activation ID quantities in the table by editing the value in the Total column if needed.

8 When all Activation ID's and quantities are finalized, click **Generate Activation Request** to create an Activation Request file.

Entitlement ID	Activation ID	Product Name	Subscription Expiration (yyyy-MM-dd)	Total
	ad1b-			1

Generate Activation Request

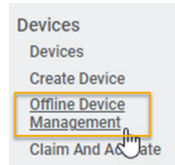
The OfflineActivationRequestX.bin file is added to the browser's default Downloads folder.



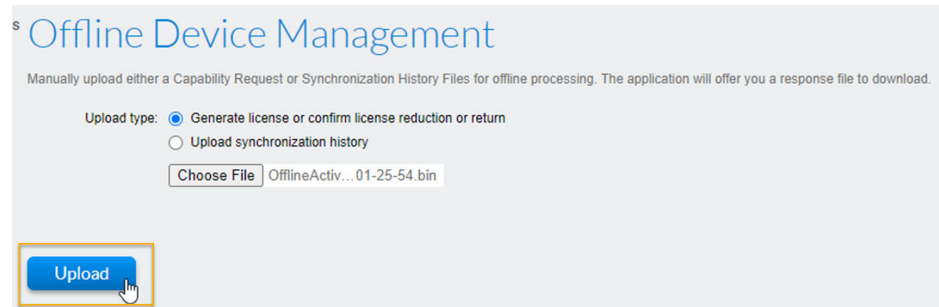
Installation

Offline Activation

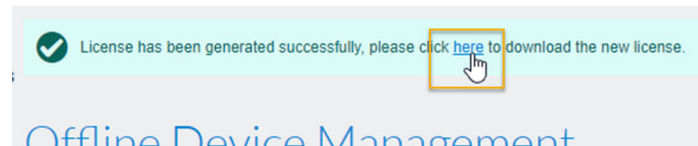
- 9 Transfer the *OfflineActivationRequestXXX.bin* file generated to a PC with Internet access.
- 10 Log into the License and Delivery Portal using your Agilent.com credentials.
- 11 Select **Offline Device Management**.



- 12 Click **Choose File**, browse to the folder where the *OfflineActivationRequest.bin* file is located and select the file.
- 13 Click **Upload**. The license is generated for download.



- 14 Click the **here** link to download the license file, *capabilityResponse.bin*, to the default download folder.



- 15 Copy the *capabilityResponse.bin* file and move it to the Installation PC.

Installation

Offline Activation

16 In the Agilent Licensing window, select the **Install Activation Response** tab.

Generate Activation Request **Install Activation Response** Activation Confirmation

Step 2 Install Activation Response file:

- Transfer and upload this *Activation Request* file to the Agilent License Portal in the "Offline Device Management" section, using the "Generate license" option. Then download the *Activation Response* file (capabilityResponse.bin) from the Agilent License Portal.
- Transfer the *Activation Response* file back to this device and browse for it. Click on the **Install Activation Response** button to upload the *Activation Response* file to this device.

Browse for the Activation Response file. *

capabilityResponse.bin

Install Activation Response

17 Browse for the Activation Response File and click **Open**.

Click **Install Activation Response**.

The software is now activated.



3

Getting Started with MassHunter Explorer Software

Now that you have the software downloaded and activated, it's time to learn more about the MassHunter Explorer software.

Next Steps

User Documentation



Data analysis and library management documentation can be found by scanning the code or navigating to <https://aglt.co/DALibMgmtDocs>. To get started, download the *MassHunter Explorer 1.0 Introduction Workbook*. View the introduction videos and use the workbook as a companion.

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>

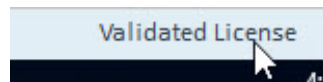
4

Transferring MassHunter Licenses

This chapter describes the preparation steps to transfer a license for MassHunter Explorer software.

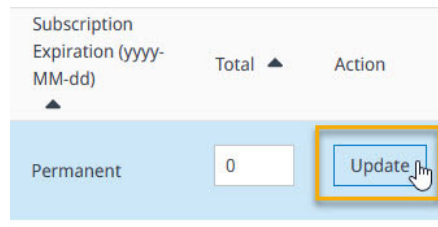
Returning MassHunter Explorer software licenses and uninstalling

- 1 Open the MassHunter Explorer software and click **Validated License** in the lower right corner of the window.



The Agilent Licensing Portal opens in a browser window.

- 2 Locate the license for the installation in the table, change the Total value to 0, then click **Update**.



The following message will flash on the screen.

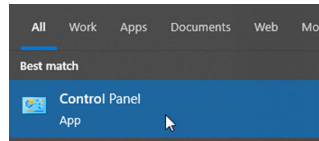


- 3 Close the MassHunter Explorer software application.

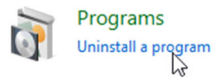
Transferring MassHunter Licenses

Returning MassHunter Explorer software licenses and uninstalling

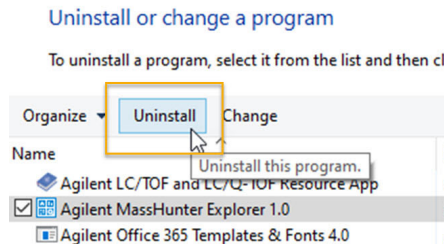
- 4 Open the Microsoft Control Panel program.



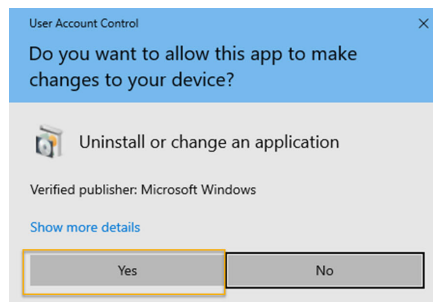
- 5 Click **Uninstall a program**.



- 6 Find and click to select **Agilent MassHunter Explorer 1.0**. The version number is listed after the program.
- 7 Click **Uninstall**.



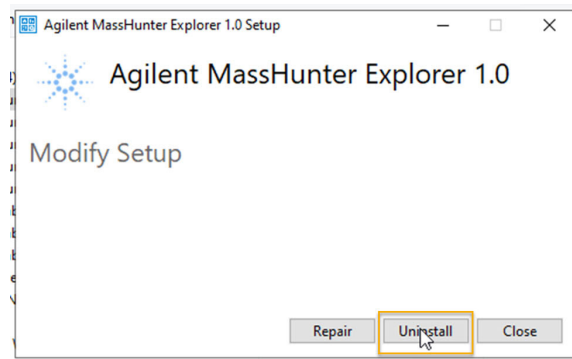
- 8 Click **Yes** to Do you want to allow this app to make changes to your device?



Transferring MassHunter Licenses

Returning MassHunter Explorer software licenses and uninstalling

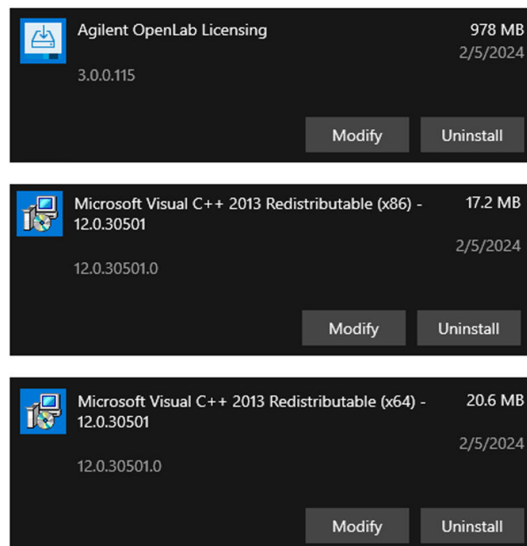
- 9 Click **Uninstall** to modify the setup.



- 10 Once the uninstallation is complete, reboot the computer.

NOTE

The MassHunter Explorer software's dependencies (.NET Framework and the License Server) are not be uninstalled with this process. If these items also need to be uninstalled (not recommended) then they can be uninstalled through the 'Add/Remove Apps' workflow available with Windows.



www.agilent.com

© Agilent Technologies, Inc. 2024
D0109800
February 2024 A.01

