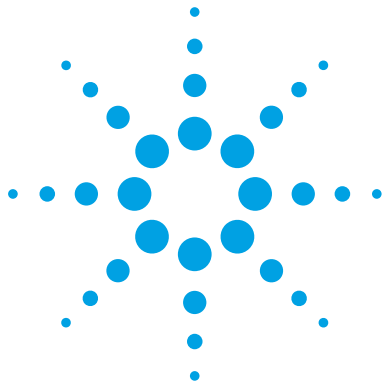




Agilent MassHunter Workstation – Offline Qualitative and Quantitative Analysis for LC/MS

Installation Guide

- Step 1. Prepare for installation 2
- Step 2. Install Microsoft Excel software 3
- Step 3. Install the DA Reprocessor - Offline Utilities 5
- Step 4. Install the Qualitative Analysis program 6
- Step 5. Install the Quantitative Analysis program 8
- Step 6. Install Quantitative Compliance program 12
- Step 7. Install Quantitative Analysis Reporting 13
- Step 8. Install Quantitative Analysis Supplemental files 16
- Step 9. Configure Excel for MassHunter 17

Use this guide to install or upgrade MassHunter Workstation software.

This guide applies to and supports only:

- MassHunter Qualitative Analysis B.07.00 and later, until superseded
- MassHunter Quantitative Analysis B.07.00 and later, until superseded
- 64-bit Windows 7 Professional SP1
- Excel 2013 and Office 2013

Do not install on Windows XP or Windows Vista.

CAUTION

MassHunter and Microsoft Excel software must be installed and set up in the order described in this guide, or the printing of reports can fail.



Agilent Technologies

Step 1. Prepare for installation

CAUTION

Do not remove programs that need only service pack installation. Remove only the programs that you plan to upgrade at this time.

- 1 Use the **Uninstall a program** or **Programs and Features** command in Control Panel to remove programs that you plan to upgrade, if they are installed on your system.

- Microsoft Excel, if you need to upgrade. Only Excel 2013, either as a standalone program or as part of the Office 2013 suite, is supported.
- Agilent MassHunter Offline Utilities
- Agilent MassHunter Workstation Qualitative Analysis
- Agilent MassHunter Workstation Quantitative Analysis
- Agilent MassHunter Workstation Quantitative Analysis Compliance
- Agilent MassHunter Workstation Quantitative Analysis Reporting

Some of these programs require a reboot when you remove them. You can wait until all of the programs are removed before you reboot your computer.

- 2 Run Windows Update to make sure you have the latest critical updates and security fixes.

- Click **Start > Control Panel > System and Security > Windows Update**.

Verify Windows Update is completed before you continue.

CAUTION

If you do not run Windows Update, your system may be vulnerable to security problems. Confirm that the LAN power management has not been started by a Windows Update program.

Step 2. Install Microsoft Excel software

Do this step if you want to upgrade and configure Microsoft Excel software to run with your MassHunter program.

MassHunter supports only Microsoft Excel 2013, or Microsoft Office 2013.

- 1** Close all MassHunter Workstation programs.
- 2** Make sure that the DA Reprocessor program is not running.
- 3** Start the installation:
 - a** Follow the instructions that came with the Microsoft Excel or Microsoft Office software or Product Key to start the installation.
 - b** If you are prompted to enter the Product Key, do so.

If you do not enter the Product Key now, Microsoft Excel reminds you later, which can interfere with report automation.

The Product Key must be verified to activate your Microsoft product. Automatic verification requires an Internet connection. If your system does not have access to the Internet, the Activation Wizard starts and allows you to verify by telephone.

- c** Click **Customize**.

Step 2. Install Microsoft Excel software

- 4 Make sure that these options, which are required by MassHunter to run reports, are enabled for installation. If not, click the program name and click **Run all from My Computer**.
 - Microsoft Office Excel
 - Office Shared Features > Visual Basic for Applications
 - Office Tools

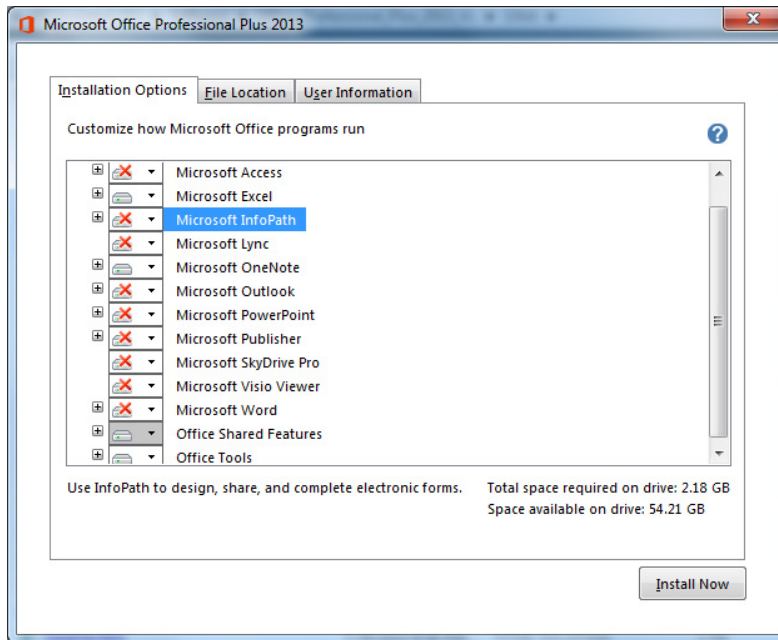


Figure 1 Microsoft Office installation options. For Microsoft Excel, you see only Microsoft Office Excel, Office Shared Features and Office Tools.

- 5 Click **Install Now** and continue the installation as instructed.
- 6 If you are prompted to get updates, click **Go to Office Online**. Follow the instructions to update Microsoft Excel or Microsoft Office.

Step 3. Install the DA Reprocessor - Offline Utilities

Do this step only if you want to reprocess LC/MS data on an offline MassHunter system.

The Offline Utilities disc contains the DA Reprocessor program. The DA Reprocessor program lets you reprocess one or more data files by running a worklist which starts the Qualitative Analysis program for each file and method in the worklist.

- 1** Insert the MassHunter Acquisition Offline Utilities disc into the disc drive.
- 2** Double-click **Setup.msi**.
- 3** In the **License Agreement** page, click **I accept the terms of the license agreement**, and then click **Next**.
- 4** In the **Choose Destination Location** page, click **Next** to accept the default destination location.

The default is **C:\Program Files (x86)\Agilent\MassHunter\Workstation\Acquisition Offline**.

- 5** In the **Installation Complete** message, click **Finish**.

Refer to the Data Acquisition online Help for more information on running the DA Reprocessor program.

Step 4. Install the Qualitative Analysis program

Do this step if you want to install the Qualitative Analysis program. If you are installing just the service pack, go directly to [step 13](#).

- 1 Verify that no other program is running on your system, including Windows Update.

Installation of the Qualitative Analysis program can fail if Windows Update is running, even in the background.

- 2 Insert the Qualitative Analysis installation disc into the DVD drive.
- 3 From the installation disc, right-click **setup.exe**, and then click **Run as administrator**.
- 4 If **Windows User Account Control** asks you to allow the program to make changes to your system, click **Yes**.
- 5 In the **Welcome** page, click **Next**.
- 6 In the **License Agreement** page, click **I accept the terms of the license agreement**, and then click **Next**.
- 7 In the **Choose destination location** page, click **Next** to accept the default destination location.

The default is **C:\Program Files\Agilent\MassHunter\Workstation\Qual**.

- 8 In the **Choose location for customer data files** page, click **Next** to accept the default destination location.

The default is **D:\MassHunter**. Click **Change** if you need to change the location.

Note that some default worklists and methods expect customer data to be in **D:\MassHunter**. If the customer data folder is set to a different location, users can get an error indication when they open these methods and worklists.

- 9 In the **Ready to install** page, review the settings, and then click **Install**.
- 10 If you get a message that indicates that the screen resolution is not set to the required resolution, click **OK**.
- 11 If you see a Notepad window that prompts you to install other options, close the Notepad window.
- 12 In the installation completion message, click **Finish**.

Step 4. Install the Qualitative Analysis program

13 If the folder **Service Packs** exists on the installation disc, then:

- 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**.

Follow the instructions to install the service pack.

Step 5. Install the Quantitative Analysis program

Do this step if you want to install the Quantitative Analysis program. If you are installing just the service pack, go directly to [step 17](#).

If Quantitative Analysis B.07.00 is already installed and you want to install only the Compliance program, continue at [“Install Quantitative Compliance program”](#) on page 12.

- 1 Verify that no other program is running on your system. Installation of the Quantitative Analysis program fails if Windows Update is running, even in the background.
- 2 Insert Quantitative Analysis Disk 1/2 into the DVD drive.
- 3 From the installation disc, right-click **setup_quant.exe**, and then click **Run as administrator**.
- 4 If **Windows User Account Control** asks you to allow the program to make changes to your system, click **Yes**.
- 5 When prompted, select the language for the installation.
- 6 If prompted, select **Install** in the Agilent Installation Qualification Tool dialog box.
- 7 In the **Welcome** page, click **Next** to continue with the installation.
- 8 In the **License Agreement** page, click **I accept the terms in the license agreement**, and then click **Next**.
- 9 In the **Destination Folder** page, click **Next** to accept the default destination location.

The default is **C:\Program Files\Agilent\MassHunter\Workstation\Quant**.

- 10 In the **Customer Data Folder** page, click **Next** to accept the default destination location.

The default is **D:\MassHunter**. Click **Change** if you need to change the location.

For a typical installation, use the same data file location that is used by the MassHunter Data Acquisition Workstation to store acquired data.

Note that some default worklists and methods expect customer data to be in **D:\MassHunter**. If the customer data folder is set to a different location, users can get an error indication when they open these methods and worklists.

11 In the **Setup Type** page, click **Typical**, **Complete**, or **Custom** setup, and then click **Next**.

- A **Typical** installation installs the basic Quantitative Analysis software.
- A **Complete** installation installs the basic Quantitative Analysis software, Quantitative Analysis Compliance software and the TPH and Blank Subtraction add-ins.
- A **Custom** installation presents you with an additional Custom Setup page that allows you to view the program components and choose which to install.

12 In the **Ready to Install the Program** page, click **Install** to begin.

When the installation begins, the **Installing Agilent MassHunter Workstation Quantitative Analysis** page appears.

13 In the **Choose Icons** page, mark the check boxes for the instruments that apply. Clear the check boxes for the instruments that do not apply. See [Figure 2](#).

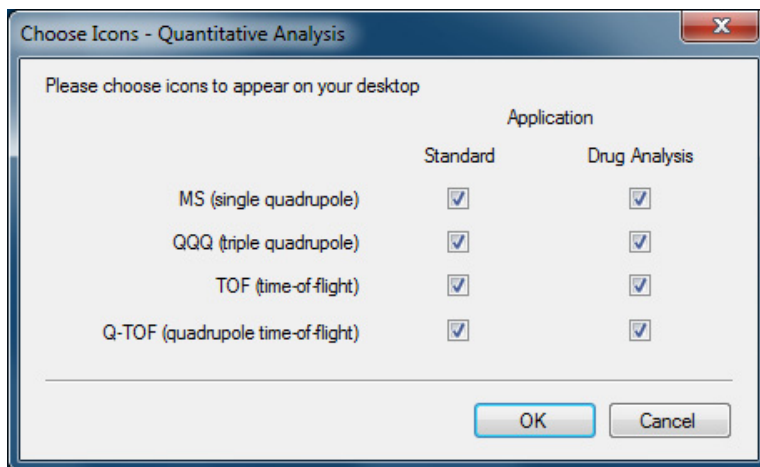


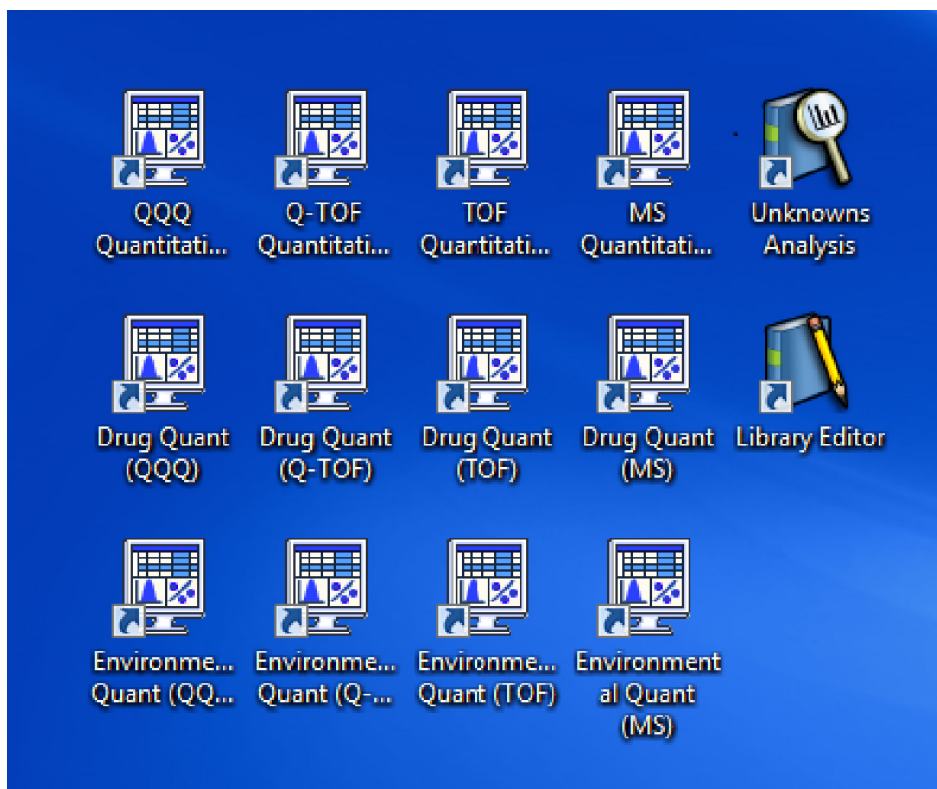
Figure 2 Choose Icons dialog box

After installation, you can change which icons are displayed. Click **Start > All Programs > Agilent > MassHunter Workstation > Quant Tools > Setup Desktop Icons**.

14 Click **OK**.

15 In the **Completed** message, click **Finish**.

16 Confirm that the icons you selected, as well as **Library Editor** and **Unknowns Analysis**, appear on your desktop.



17 If a service pack is available:

- a** If the folder **Service Packs** exists on the installation disc, then open the **Service Packs** folder. If the folder contains another folder, such as **SP1**, open that folder.

Step 5. Install the Quantitative Analysis program

As an alternative, the service pack file can be provided to you by an Agilent representative or support service.

- b** Locate the installation file, ending with **SPx.exe**.
- c** Right-click the installation file and click **Run as administrator**.
- d** Follow the instructions to install the service pack.

Step 6. Install Quantitative Compliance program

Do this step if you need to enable compliance for the Quantitative Analysis program.

- 1** From the Quantitative Analysis installation disc, right-click **setup_quantcompliance.exe** and click **Run as administrator**.
- 2** Select desired language and click **OK**.
- 3** In the installation **Welcome** page, click **Next** to continue.
- 4** In the License Agreement section, click **I accept the terms of the license agreement**, and then click **Next**.
- 5** Click **Install**.
- 6** Click **Finish** when install is complete.

Step 7. Install Quantitative Analysis Reporting

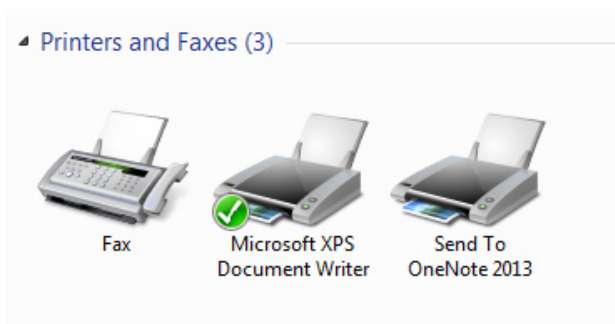
Do this step only if you need to upgrade your Quantitative Analysis Reporting program. If you are installing just the service pack, go directly to [step 10](#).

- 1 Verify that no other program is running on your system. Installation of the Quantitative Analysis Reporting program will fail if Windows Update is running, even in the background.
- 2 From the root directory of the Quantitative Analysis installation disc, right-click **setup_quantreport.exe**, and then click **Run as administrator**.
- 3 Select desired language from the drop-down menu and click **OK**.
- 4 In the **Welcome** page, click **Next** to continue with the installation.
- 5 In the **License Agreement** page, mark the **I accept the terms of the license agreement**, and then click **Next**.
- 6 If needed, in the **Destination Folder** page, click **Change** to change the default destination location for the report templates. The default is **D:\MassHunter\Report Templates\Quant**.
- 7 In the **Setup Type** page, click **Typical**, **Complete**, or **Custom** setup, and then click **Next**.
 - A **Typical** or **Complete** installation installs all the program components and requires the most disk space.
 - A **Custom** installation presents you with an additional Custom Setup page that allows you to view the program components and choose which to install.
- 8 In the **Ready to Install the Program** page, click **Install** to begin.

When the installation begins, the **Installing Agilent MassHunter Workstation Quantitative Analysis Reporting** page appears.
- 9 In the **Completed** page, click **Finish**.
- 10 If a service pack is available:
 - a If the folder **Service Packs** exists on the installation disc, then open the **Service Packs** folder. If the folder contains another folder, such as **SP1**, open that folder.

As an alternative, the service pack file can be provided to you by an Agilent representative or support service.

- b** Locate the installation file, ending with **Reporting_SPx.exe**.
 - c** Right-click the installation file and click **Run as administrator**.
 - d** Follow the instructions to install the service pack.
- 11** Make sure the appropriate print drivers are installed.
- To use the HP LaserJet P3005d or P3015d printer, install the print driver from <http://welcome.hp.com/country/us/en/support.html>.
 - To print reports in the A4 format to HP PCL6 printers, install the printer driver patch from <http://support.microsoft.com/kb/935843>.
 - To use the Canon LBP3310 printer for Japanese Windows, install the printer driver from the installation disc.
- 12** Set the Microsoft XPS Document Writer as the default printer:
- a** Click **Start > Devices and Printers**.
 - b** Right-click **Microsoft XPS Document Writer** and click **Set as Default Printer**.



CAUTION

If no printer is set up as the default printer, reporting can fail.

If you installed Excel 2013 or Office 2013, you must change the default printer to **Microsoft XPS Document Writer**. These products automatically install Microsoft OneNote and set the default printer to OneNote. OneNote breaks some Excel report formats and causes extra pages to be printed.

Step 7. Install Quantitative Analysis Reporting

The Microsoft XPS Document Writer formats the Excel workbooks three to four times faster for networked printers because the formatting page set-up information is stored on the local computer. The report is processed by the Microsoft XPS Document Writer, but it is actually printed on the printer selected in the MassHunter program.

Step 8. Install Quantitative Analysis Supplemental files

Do this step if you want to copy these to your computer from the installation disc:

- Manuals
- Report template index
- Technical videos outline
- Data

- 1** From the MassHunter Quantitative Analysis installation disc, open the **Supplemental** folder. Right-click **Setup_Supplemental.exe**, and then click **Run as administrator**.
- 2** If **Windows User Account Control** asks you to allow the program to make changes to your system, click **Yes**.
- 3** Select desired language from the drop-down menu and click **OK**.
- 4** In the **Welcome** page, click **Next** to continue with the installation.
- 5** In the **License Agreement** page, mark the check box for **I accept the terms of the license agreement**, and then click **Next**.
- 6** If needed, in the **Destination Folder** page, click **Change** to change the default destination location for the supplemental items. The default is **D:\MassHunter**. Click **Next**.
- 7** In the **Setup Type** page, click **Complete** or **Custom** setup, and then click **Next**.
 - A **Complete** installation installs all the available components and requires the most disk space.
 - A **Custom** installation presents you with an additional Custom Setup page that allows you to view the program components and choose which to install.
- 8** In the **Ready to Install the Program** page, click **Install** to begin.
- 9** In the **Completed** page, click **Finish**.
- 10** To view the supplemental materials, click **Start > All Programs > Agilent MassHunter Workstation > Quant Supplemental**.

Step 9. Configure Excel for MassHunter

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

- 1** Click **Start > Control Panel > Programs and Features**.
- 2** If you need to configure Excel for Qualitative Analysis:
 - a** Right-click **Agilent MassHunter Workstation Qualitative Analysis**.
 - b** Click **Change**, and then click **Repair**.
 - c** Follow the instructions to complete the process.
- 3** If you need to configure Excel for Quantitative Analysis:
 - a** Right-click **Agilent MassHunter Workstation Quantitative Analysis**.
 - b** Click **Change**, and then click **Repair**.
 - c** Follow the instructions to complete the process.

www.agilent.com

In This Book

This book contains installation instructions to install or upgrade an offline copy of the Agilent MassHunter Workstation Software.

Microsoft®, Windows 7®, Windows XP®, Windows Vista®, and Excel® are registered trademarks of Microsoft Corporation in the US and/or other countries.

© Agilent Technologies, Inc. 2014

Printed in USA
Revision A, December 2014



G3335-90181



Agilent Technologies