

Mass Profiler Professional

Installation Guide

- To do a new installation 2
- To update from MPP 13.0 2
- To update from MPP 13.1 3
- To update from MPP 13.1.1 3

Before you install Mass Profiler Professional (MPP) software, make sure that compatible software for your application is installed.

Compatibility

Table 1 Software compatibility

Application	Compatible software
LC/MS data extraction	<ul style="list-style-type: none">MassHunter Profinder B.06.00 SP1 or higher.MassHunter Qualitative Analysis B.05.00 or later with 64-bit Windows 7.
GC/MS data extraction (MassHunter)	<ul style="list-style-type: none">MassHunter Qualitative Analysis B.05.00 or later with 64-bit Windows 7.Data must first be converted from the MSD ChemStation data format into the MassHunter data file format via the translator which is provided with the MassHunter for GC/MS products
GC/MS data extraction (ChemStation)	<ul style="list-style-type: none">G1701EA MSD ChemStation E.02.00 SP2 or higher
Feature finding in GC/MS data acquired in the MSD ChemStation format	<ul style="list-style-type: none">NIST AMDIS software required



Table 1 Software compatibility (continued)

Application	Compatible software
Identification of GC/MS features via EI-MS library search	<ul style="list-style-type: none"> G1033A NIST 11 software can optionally be purchased and installed. Older revisions of the NIST library will also work.
Identification of endogenous metabolites in GC/MS data files via EI-MS library search	<ul style="list-style-type: none"> G1676AA Fiehn GCMS Metabolomics RTL Database software can optionally be purchased and installed.
Identification of endogenous metabolites in LC/MS data files via accurate mass database search	<ul style="list-style-type: none"> Agilent G6825AA Personal METLIN Metabolite Database or G3874AA MassHunter METLIN Metabolite PCDL software can optionally be purchased and installed. Version B.05.00 is recommended. MassHunter PCDL Manager B.07.00 SP1 or higher is required.

To do a new installation

- From the installation media, run **MPP14.0_Setup.exe** as administrator. To do so, right-click **MPP14.0_Setup.exe** and select **Run as administrator**.

If you do not have a valid Order ID, you can get a temporary license at <http://lcosgens.cos.agilent.com/gsLicense/MPPTempLicense.html>. Refer to the *Quick Start Guide* for more information about license activation for Mass Profiler Professional and Pathway Architect.

To update from MPP 13.0

- Update to MPP 13.1.1:
 - Start the Mass Profiler Professional 13.0 program.
 - Select **Help > Update Product > From Agilent Server**.
If you cannot update from the Agilent server, click **Help > Update Product > From file**. From the Supplemental disc, in the **Updates** folder, select **MPP13.1.update**.
- Continue at “[To update from MPP 13.1.1](#)”.

To update from MPP 13.1

1 Update to MPP 13.1.1:

- a** Start the Mass Profiler Professional 13.1 program.
- b** Select **Help > Update Product > From Agilent Server**.

If you cannot update from the Agilent server, click **Help > Update Product > From file**. From the Supplemental disc, in the **Updates** folder, select **MPP13.1.1.update**.

2 Continue at “[To update from MPP 13.1.1](#)”.

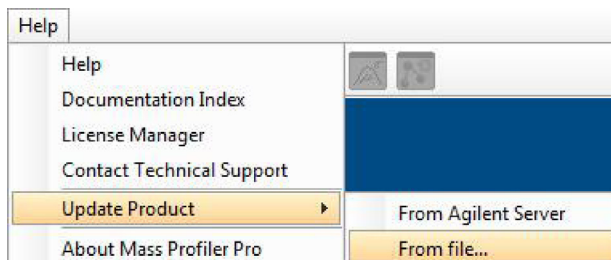
To update from MPP 13.1.1

This update removes MPP 13.1.1 from your system.

Do not use any other program or tool to remove MPP 13.1.1 before you install the MPP 14.0 update. You must update from the update file on the MPP 14.0 Supplemental disc.

Note that the MPP 14.0 update is available only to customers who have purchased an upgrade product or have an active Software Maintenance Agreement (SMA).

- 1** Start the Mass Profiler Professional 13.1.1 program.
- 2** Select **Help > Update Product > From file**.
- 3** From the Supplemental disc, in the **Updates** folder, select **MPP14.0.update**.



The conversion of projects to the new MPP 14.0 format can take 30 minutes to an hour, depending upon the number of projects.

www.agilent.com

In This Guide

This guide contains instructions to install or upgrade to the Mass Profiler Professional 14.0 software.

Windows 7® is a registered trademark of Microsoft Corporation in the US and/or other countries.

This guide applies to Mass Profiler Professional 14.0.

© Agilent Technologies, Inc. 2015

Revision A, November 2015



G3835-90024



Agilent Technologies