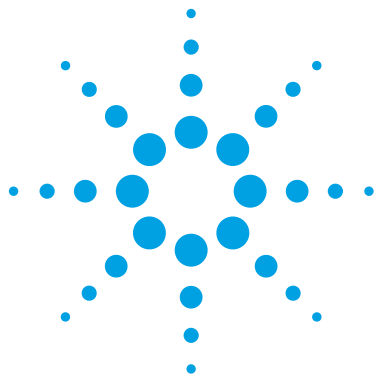




Agilent G3835AA MassHunter Mass Profiler Professional Software

Quick Start Guide

1. Getting Started 6
2. User Interface 10
3. Basic Tasks 15
4. Advanced Tasks 24
5. Acknowledgment and Citations 24

What is Agilent Mass Profiler Professional?

Agilent Mass Profiler Professional (MPP) software is a powerful chemometrics platform designed to exploit the high information content of mass spectra (MS) data and can be used in any MS-based differential analysis to determine relationships among two or more sample groups and variables. MPP provides advanced statistical analysis and visualization tools for GC/MS, LC/MS, CE/MS, and ICP-MS data analysis. MPP also integrates smoothly with Agilent MassHunter Workstation, Spectrum Mill, and ChemStation software and is the only platform that provides integrated identification/annotation of compounds and integrated pathway analysis for metabolomic and proteomic studies. The system also enables automated sample class prediction that revolutionizes mass spectrometer-based qualitative analysis of unknown samples in many applications. MPP is ideally suited for applications characterized by complex sample matrices such as metabolomics, proteomics, natural products, food, beverages, flavors, fragrances, and environmental analyses.



Agilent Technologies

Features of Mass Profiler Professional

- Allows differential analysis of two or more sample sets from one or multiple MS analysis platforms in a single project.
- Designed for both mass spectrometrists and statisticians, with guided and advanced workflows.
- Provides comprehensive statistical tools including ANOVA, PCA, volcano plots, hierarchical trees, SOMs, QT clustering, and five different methods for class prediction.
- Generates inclusion lists for subsequent MS/MS-based analyses to facilitate identification of compounds of interest.
- The ID Browser allows integrated identification and annotation of compounds using LC/MS Personal Compound Database Libraries and GC/MS libraries (NIST and Agilent Fiehn Metabolomics).
- Allows import and export of data in several formats e.g. to further extend and customize statistical analysis and visualization capabilities using R-Scripts (see *Agilent GeneSpring GX R-integration package - Installation and Application Guide*).
- Integrates smoothly with Spectrum Mill to enable TOF/Q-TOF based discovery and QQQ-based verification of protein biomarkers.
- Integration with GeneSpring allows the correlation of pathways between MS (metabolomics, proteomics) and microarray data (transcriptomics) in one project, enabling true integrated biology studies.
- Class prediction models from GC/MS, LC/MS, CE/MS and ICP/MS data can be used with the new, standalone Automated Sample Class Prediction software to enable high-throughput routine classification of unknown samples.

What's new in MPP 14.0?

- New analytical capabilities
 - Analysis of *Agilent Proteomics Results* (APR) files exported from SpectrumMill and Skyline Software. The tool supports results for both label free and labeled experiments (SILAC, iTRAQ, TMT).
 - Quilt Plot views, miniature heatmap data overlays, are available when viewing pathways to provide enhanced experiment data visualization.
 - Statistical analysis on proteins, peptides, and post-translational modifications of proteins and peptides.
 - Support for stable isotope label experiments (SILAC, ICAT, iTRAQ, TMT) via Spectrum Mill.
 - Peptide and Protein level analyses and filters in:
 - Quality Control - Filter on Peptides and Filter on Proteins
 - Spreadsheet views - sort by peptide sequence
 - Workflows - Filter by Peptide Profile and Filter by Protein Modifications
 - Statistical and correlation analyses include peptide and protein results
 - Visualize peptide and protein results in various plots
 - Visualize peptide and protein results on pathways
 - Export proteins, genes, and metabolites from selected pathways.
 - Annotated experimental results on KEGG Pathway maps.
- Usability enhancements
 - Legend for pathways is added on the side panel and is now visible during full view mode.
 - Duplicate pathways, created during project import, do not appear during SEA and MOA pathway analyses.
 - Ability to enable and disable automatic check for product and technology updates during software startup.
 - Multiple colors are used to denote the experiment type for heatstrips drawn on pathway views.
 - Dendrogram view is updated to respond dynamically to changes in the active entity list.
 - Ability to export PCA plot images in SVG and EPS formats.

- Export of entities in cluster order.
- BioPAX file import allows special characters in the pathway name.
- Imported BioPAX pathways as BioCyc pathways are available for analysis.
- MassHunter complementary features
 - MassHunter ID Browser B.07.00 Service Pack 1

Where to Find More Information

Online help

Press F1 To get more information about a pane, window, or dialog box place the cursor on the part of the pane, window, or dialog box of interest and press **F1**.

Help Menu Click **Help > Documentation Index** to access the release notes, quick start guides, and HTML and PDF versions of the *Mass Profiler Professional User Manual*.

Documents

Mass Profiler Professional

- Mass Profiler Professional Familiarization Guide (p/n G3835-90026)
- Mass Profiler Professional Application Guide (p/n G3835-90028)
- Agilent MassHunter Mass Profiler Professional User Manual.

You can find a PDF copy in the MPP installation folder

C:\Program Files\Agilent\MassHunter\Workstation\Mass Profiler Professional\docs\manual\MS, if you installed MPP using the default directory structure.

Workflow Guides and Overviews

- Agilent Metabolomics Workflow - Discovery Workflow Guide (p/n 5990-7067EN)
- Agilent Metabolomics Workflow - Discovery Workflow Overview (p/n 5990-7068EN)
- Integrated Biology with Agilent Mass Profiler Professional - Workflow Guide (p/n 5991-1909EN)
- Integrated Biology with Agilent Mass Profiler Professional - Workflow Guide Overview (p/n 5991-1910EN)
- Class Prediction with Agilent Mass Profiler Professional Workflow Guide (p/n 5991-1911EN)
- Class Prediction with Agilent Mass Profiler Professional Workflow Overview (p/n 5991-1912EN)

MassHunter Qualitative and Quantitative Analysis

- MassHunter Qualitative Analysis Familiarization Guide for LC/MS
- MassHunter Qualitative Analysis Familiarization Guide for GC/MS
- MassHunter Quantitative Analysis Familiarization Guide

Training

Road Show Slide Presentation Metabolome Analysis - From Sample Prep Through Data Analysis (Metabolomics Road Show, March 2011)

Training Videos Training videos can be found on the Supplemental disc.

Training Courses Visit www.chem.agilent.com to view a listing of training courses for MPP.

1. Getting Started

How do I get started?

This *Quick Start Guide* helps you launch MPP, activate your license, review the MPP user interface, and create a project and an experiment that you import data into and then use to begin your analysis. If you are already familiar with the MPP software, after you install MPP you can immediately start using the preloaded demonstration project called “Malaria.”

Start and Explore Mass Profiler Professional

- 1 Double-click the **Mass Profiler Professional** icon  located on your desktop, or click **Start > All Programs > Agilent > MassHunter Workstation > Mass Profiler Professional > Mass Profiler Professional**.
- 2 Activate your license.
The first time you start MPP you are prompted to enter your **Mass Profiler Professional Order ID** before you can use the software (see [“Activating a License”](#) on page 7).
- 3 Review the following three sections in this *Quick Start Guide*:
[“2. User Interface”](#) on page 10
[“3. Basic Tasks”](#) on page 15
[“4. Advanced Tasks”](#) on page 24
- 4 Review the [Mass Profiler Professional Familiarization Guide \(p/n G3835-90026\)](#) (see [“Documents”](#) on page 4). This guide helps you quickly create a project, create an experiment, import and organize data, and create an initial differential expression using pre-installed demonstration data - *Malaria Demo*.
- 5 Review the [Mass Profiler Professional Application Guide \(p/n G3835-90028\)](#) (see [“Documents”](#) on page 4). This guide provides you with more detail about the MPP software and helps you create a project, create an experiment, import and organize data, and create an initial differential expression with your own data.
- 6 Perform advanced analyses to meet your project needs using the operations available in the Workflow Browser (see [“Workflow Browser”](#) on page 13).

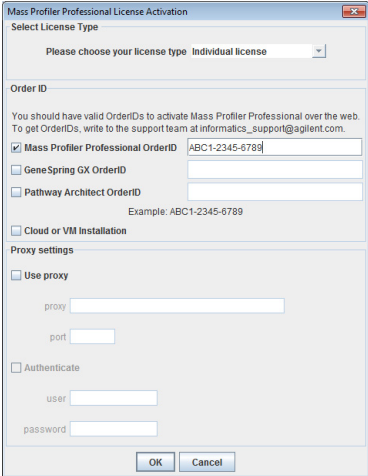
License Management

Activating a License

The first time you start MPP you enter your Order ID. License activation information is entered in the **Mass Profiler Professional License Activation** dialog box shown in [Figure 1](#). An Internet network connection is required.

- 1 Mark **Mass Profiler Professional Order ID**. If you have additional software product licenses, such as for Pathway Architect or GeneSpring GX, mark your additional products.
- 2 Type your **Mass Profiler Professional Order ID** exactly as written, for example ABC1-2345-6789. If you have additional software product licenses, such as for Pathway Architect or GeneSpring GX, type your additional product Order IDs.
- 3 Click **OK**.

The Mass Profiler Professional software automatically validates your license and starts Mass Profiler Professional.



The image shows a screenshot of the 'Mass Profiler Professional License Activation' dialog box. The window has a title bar with the text 'Mass Profiler Professional License Activation' and standard Windows window controls. The main content area is titled 'Select License Type' and contains a dropdown menu with 'Individual license' selected. Below this is the 'Order ID' section, which includes a text box containing 'ABC1-2345-6789'. There are three checkboxes: 'Mass Profiler Professional OrderID' (checked), 'GeneSpring GX OrderID', and 'Pathway Architect OrderID'. Below these is a text box for 'Example: ABC1-2345-6789'. The 'Proxy settings' section has a 'Use proxy' checkbox (unchecked) and fields for 'proxy', 'port', 'user', and 'password'. At the bottom are 'OK' and 'Cancel' buttons.

Figure 1 License Activation dialog box

1. Getting Started

Moving a License

MPP licensing is automatically enforced. You may move a license to another computer by surrendering it on the current computer and then activating it on the new computer. If you need to uninstall MPP, it is always best to surrender the existing license before un-installing MPP so that the license is available to be moved to another computer.

To surrender your license:

1 Click **Help > License Manager** in the menu bar.

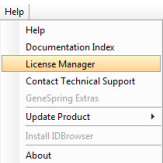


Figure 2 License Manager is available from the Help menu.

2 Click **Surrender** in the **License Manager** dialog box.

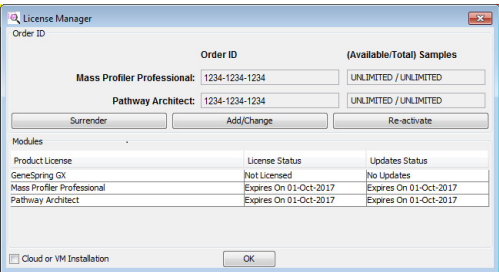


Figure 3 License Manager dialog box

3 Mark **Mass Profiler Professional** and any additional product Order IDs to surrender.

4 Click **OK**.

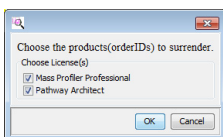


Figure 4 Order ID surrender product selection dialog box

5 Click **Yes** in the **Confirm Surrender** dialog box.

Because your license is no longer active on the current computer, MPP software immediately closes when you click **OK**.

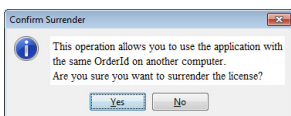


Figure 5 Confirm Surrender dialog box

6 Click **OK** in the **Surrender Status** dialog box.

Because your license is no longer active on the current computer, MPP software immediately closes.

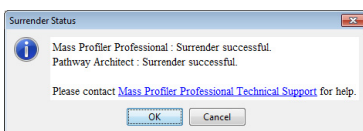


Figure 6 Surrender Status dialog box

- 7** Activate your license Order ID on the new computer (see “[Activating a License](#)” on page 7).
- 8** (Optional) Click **Start > All Programs > Agilent > MassHunter Workstation > Mass Profiler Professional > Uninstall** to remove MPP from the current computer.

2. User Interface

Main Functional Areas

The functional areas of Mass Profiler Professional are shown in [Figure 9](#) on page 11. The main MPP window consists of four parts: (1) the Menu Bar, (2) the Toolbar, (3) the Display Pane, and (4) the Status Bar.

Note: Help and detailed information regarding the various fields and statistical treatments are available by pressing **F1** or by referring to the *Mass Profiler Professional User Manual*.

1. Menu Bar The menu bar ([Figure 7](#)) provides actions that are used for managing your projects, experiments, pathways, and display pane views.

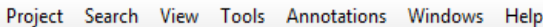


Figure 7 Menu bar

2. Toolbar The toolbar ([Figure 8](#) on page 10) is located below the menu bar and contains five sections of buttons for commonly used tasks:

Project section: New project, Open project, and Close project

Experiment section: New experiment, New MOA experiment, and Add experiment

Entity List section: Create entity list from selection, Inspect selected entity, and Import entity list (From File, From Dropbox, From Google Drive)

Plots, Statistics, and Pathways section: Scatter plot, Profile plot, Histogram plot (Single, Multiple), Matrix plot, Venn Diagram plot (Entity List, Pathway List), Box and Whisker plot, Violin plot, Heatmap plot, Data spreadsheet, Summary statistics, Create new pathway list, and Select data source for plots (Raw, Raw Scaled, Log2 Normalized, Log10 Normalized).

Sidebar section: Hide or show the sidebars that contain the Project Navigator, Workflow Browser, and Legend.

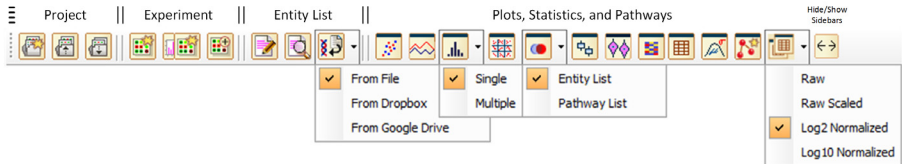


Figure 8 Toolbar and associated drop-down lists

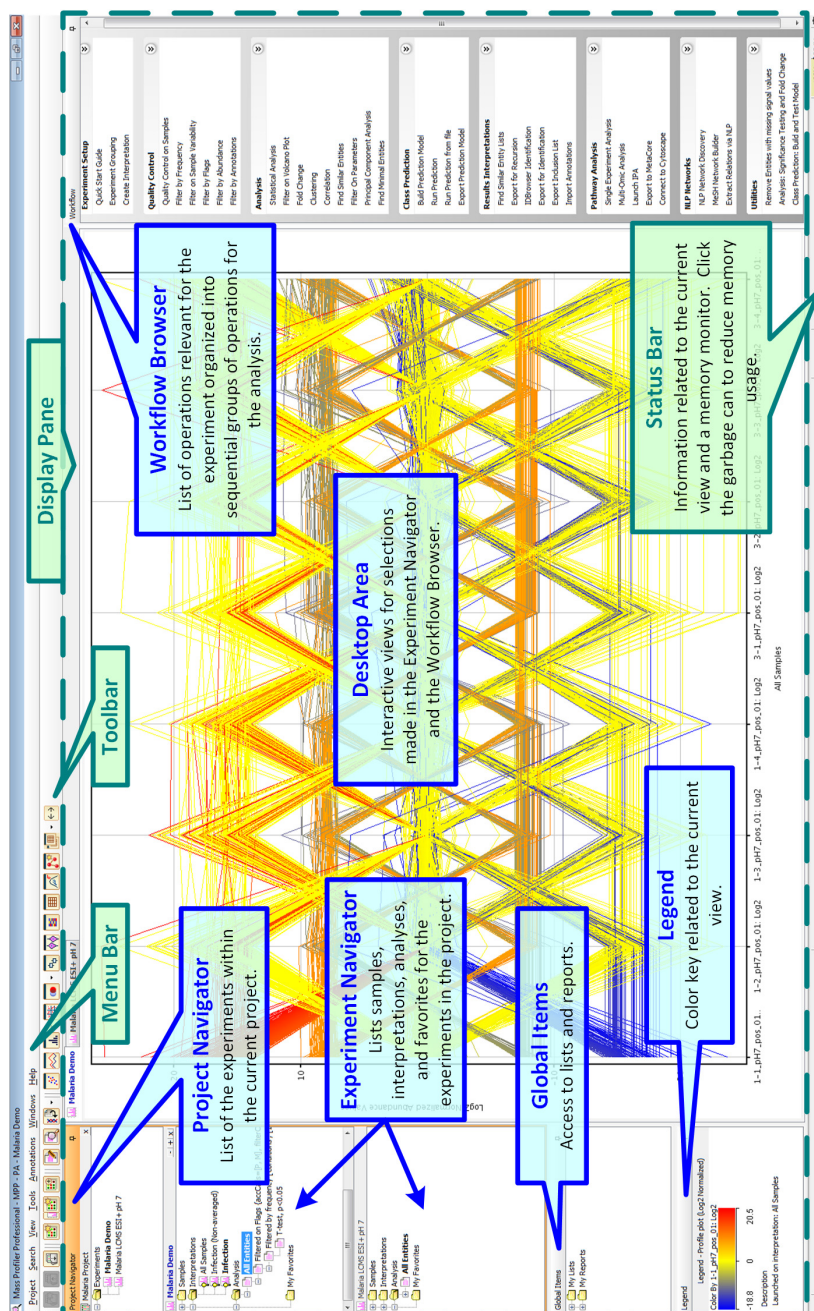


Figure 9 The main functional areas of the MPP software

Note: This figure is rotated on the page for printing.

2. User Interface

3. Display Pane The display pane, see [Figure 9](#) on page 11, is further divided into six areas – Project Navigator, Experiment Navigator, Global Items, Desktop Area, Workflow Browser, and Legend.

Project Navigator

Displays the current project and lists all the experiments with the project.

Experiment Navigator

Displays information related to Samples, Interpretations, Analysis, and My Favorites in folders related to the selected experiment in the Project Navigator. Each experiment within a project has a separate experiment navigator pane.

Global Items

Displays list and report information for your project.

Desktop Area

Display of one or more interactive views associated with the experiments. Each view in the desktop area is marked with a tab and you can configure each view separately. Click **Windows > Tile** to arrange multiple views on the desktop. The current experiment, interpretation, and analysis is noted using bold font in the Project Navigator and Experiment Navigator.

A right-click anywhere in the active view shows you a menu of options ([Figure 10](#) on page 12) to customize the view, copy the view to the system clipboard, and export the view as an image file (bmp, jpg, jpeg, png, and tiff) or as an HTML file.

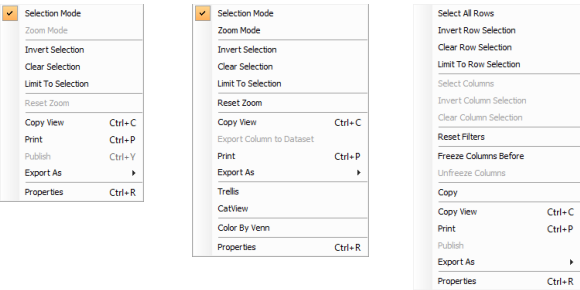


Figure 10 Right-click options within the desktop view automatically adjust to the type of view. These examples are from views containing plots and spreadsheets.

Workflow Browser

The Workflow Browser, [Figure 11](#), is organized into sequential groups of operations for the analysis of your data: Experiment Setup, Quality Control, Analysis, Class Prediction, Results Interpretations, Pathway Analysis (*optional*), NLP Networks (*optional*), and Utilities.

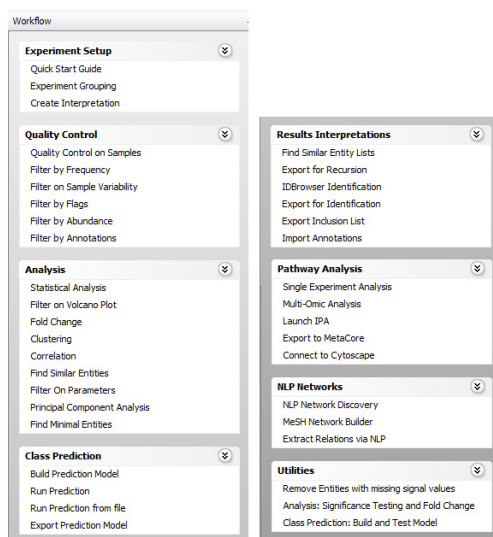


Figure 11 Workflow Browser operations

Legend

The Legend ([Figure 12](#)) shows the key (scale) to the use of color in the active desktop view. Right-click on the Legend window to copy and export the legend as an image.



Figure 12 Legend

2. User Interface

4. Status Bar The status bar (Figure 13) has three informative areas: the Status Area, the Ticker Area, and the Memory Monitor.

Status Area: Displays high-level information of the current view such as the number of rows and columns in table views and the number of entities or conditions selected in plot views.

Ticker Area: Displays the coordinates of the cursor in active plot views or the entity identification and value in table views.

Memory Monitor: Displays the total memory being used and the total memory allocated by MPP. You can click the **Garbage Can** icon  at any time to reduce the memory usage.

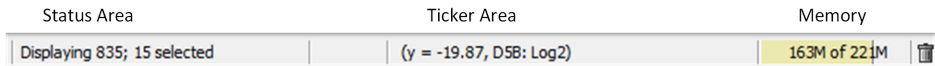


Figure 13 Status Bar

3. Basic Tasks

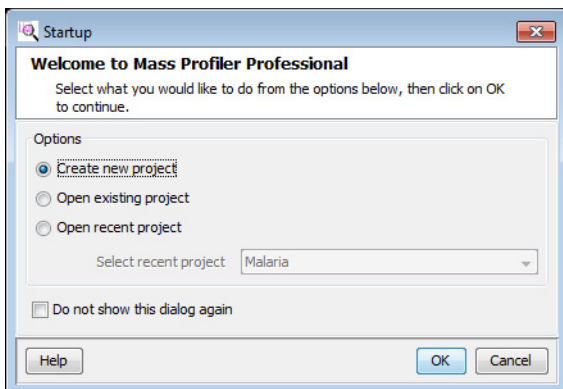
Set up a new project

A project is a “container” for a collection of experiments, each experiment contains your samples, interpretations, and analyses. A project can have multiple experiments on different sample types and organisms.

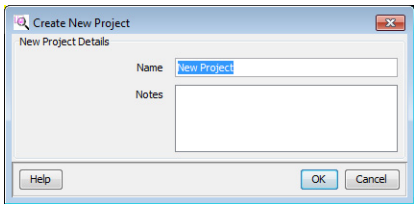
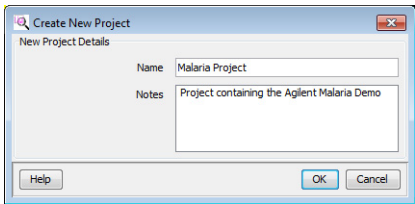
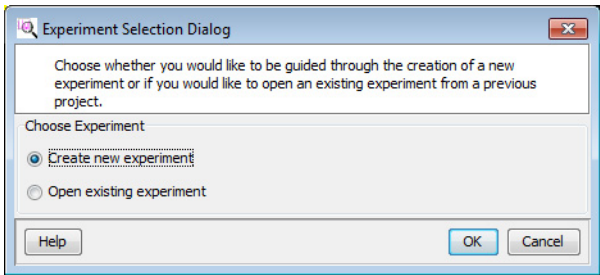
Create a new project and experiment for your data

A new project and experiment is created through four sequential dialog boxes: (1) Startup, (2) Create New Project, (3) Experiment Selection, and (4) New Experiment. The Agilent *Malaria Demo* data supplied with MPP is used as an example in each step. You may substitute the malaria information with information for your data.

Steps	Detailed Instructions	Comments
1 Start Mass Profiler Professional.	a Click the Mass Profiler Professional icon  on your desktop, or click Start > All Programs > Agilent > MassHunter Workstation > Mass Profiler Professional > Mass Profiler Professional.	The Startup dialog box helps you quickly set up a new project (this section) or continue your analysis with a previous project (see “Open a recent or existing project” on page 20).
2 Create a new project from the Startup dialog box.	a Click Create new project. b Click OK.	Create new project provides you with the option to create a new experiment or import an experiment from an existing project into the new project. - After closing an open project, you may create a new project from the Menu bar, click Project > New Project, or from the Toolbar, click the New project  button.



3. Basic Tasks

Steps	Detailed Instructions	Comments
3	Type descriptive information in the <i>New Project Details</i> section of the Create New Project dialog box.	
	a Type Malaria Project in Name. b Type Project containing the Agilent Malaria Demo in Notes. c Click OK.	• The project name and notes may be viewed and edited at any time using the Project Inspector by clicking Project > Inspect Project from the menu bar.
	 > 	
4	Create a new experiment in the Experiment Selection Dialog dialog box.	
	a Click Create new experiment. b Click OK.	• You may create or add a new experiment to your project from the Menu bar, click Project > New Experiment, or from the Toolbar, click the New experiment  button, or click the Add experiment  button. • If you click Open existing experiment, you are prompted for the experiment to add to the project as described in “Open a recent or existing project” on page 20.
		

Steps	Detailed Instructions	Comments
5	Type and select information that guides the experiment creation in the New Experiment dialog box.	
	a Type the descriptive name <i>Malaria Demo</i> in Experiment name. This name may be different from the project name. b Select Mass Profiler Professional for Analysis type. c Select Unidentified for the Experiment type. d Select Analysis: Significance Testing and Fold Change for Workflow type. e Type <i>Malaria LCMS ESI+ pH 7</i> in Experiment notes. f Click OK.	Experiment type determines how Mass Profiler Professional manages the data. Combined (Identified + Unidentified) is the proper Experiment type when you are unsure if the data has been identified in full or in part, or when other MassHunter software has been previously used to identify some of the compound features. Identified or Proteomics (Identified) is the proper Experiment type selection when the compounds have been identified by compound, formula, and/or CAS number. Unidentified is the proper Experiment type selection when the compounds have only been identified by their molecular features of neutral mass and retention time.

>

- Your new project is now set up.
- You are **immediately** guided through the steps to import and organize your data followed by the steps to create an initial analysis.

Import your data using the “MS Experiment Creation Wizard”

After you set up a project and create an experiment, the “MS Experiment Creation Wizard” guides you through the necessary steps to organize your experiment, import your data, define your experiment variables, and prepare your data for analysis. The “MS Experiment Creation Wizard” flow diagram is illustrated [Figure 14](#). The data preparation includes grouping, filtering, alignment, normalization, and baselining. The [Mass Profiler Professional Familiarization Guide \(p/n G3835-90026\)](#) (see “Documents” on page 4) guides you through importing and organizing the *Malaria Demo* data. See the [Mass Profiler Professional Application Guide \(p/n G3835-90028\)](#) for details to import non-Agilent data.

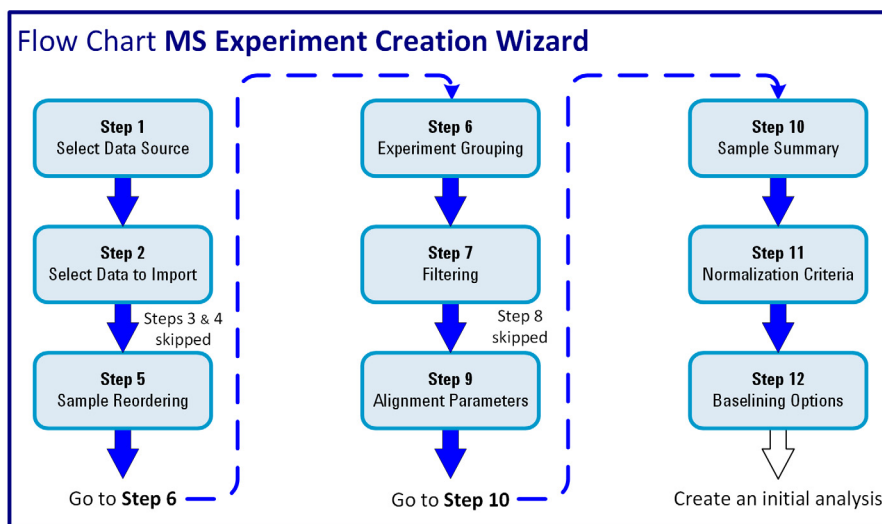


Figure 14 Steps in the “MS Experiment Creation Wizard” for the *Malaria Demo*.

Perform your initial differential analysis using the Analysis: Significance Testing and Fold Change workflow

The “Significance Testing and Fold Change Wizard” guides you through the steps necessary to enter parameters and values that improve the quality of your results and produce an initial differential expression for your analysis. The “Significance Testing and Fold Change Wizard” flow diagram is illustrated in [Figure 15](#) on page 19. The [Mass Profiler Professional Familiarization Guide \(p/n G3835-90026\)](#) (see “Documents” on page 4) guides you through an initial differential analysis using the *Malaria Demo* data.

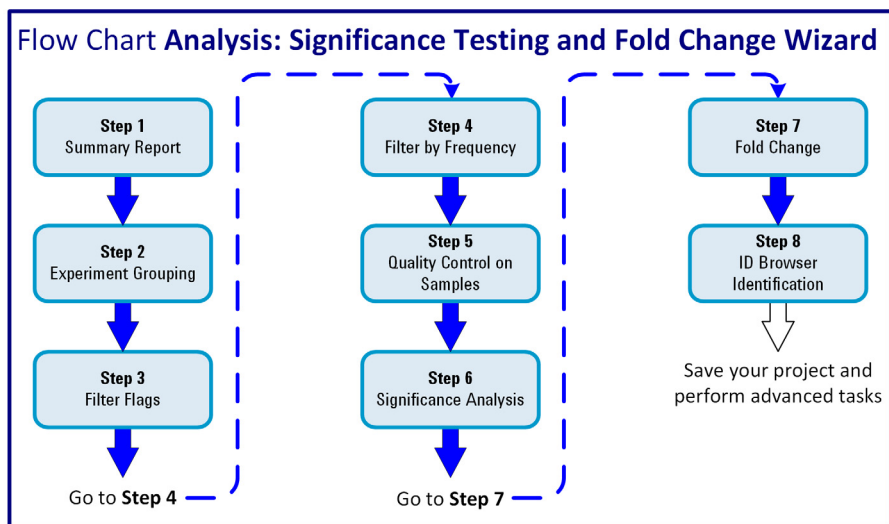


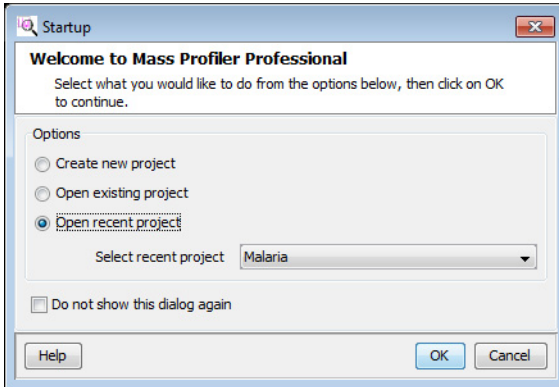
Figure 15 Steps in the "Analysis: Significance Testing and Fold Change Wizard."

Open a recent or existing project

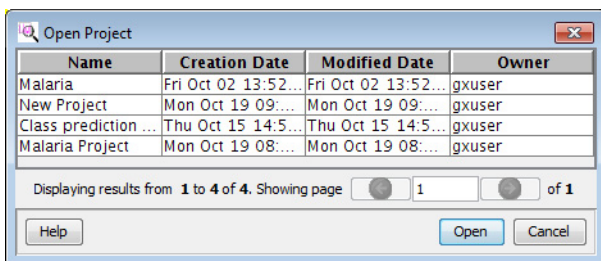
When you start Mass Profiler Professional the **Startup** dialog box allows you to quickly open a recent project, an existing project on your computer, or import an experiment from an existing project into a new project.

The steps presented are recommended if you are already familiar with the operations available in the Workflow Browser. If you are new to using Mass Profiler Professional start with “Set up a new project” on page 15 and then refer to the [Mass Profiler Professional Familiarization Guide \(p/n G3835-90026\)](#) (see “Documents” on page 4).

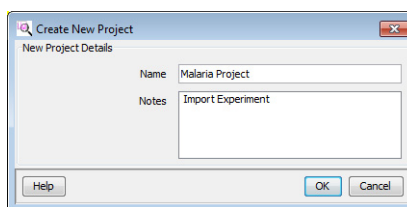
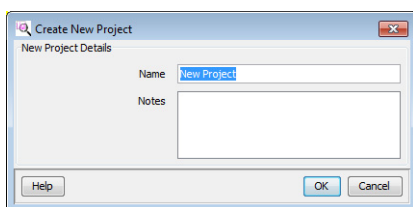
Steps	Detailed Instructions	Comments
1 Start Mass Profiler Professional.	a Click the Mass Profiler Professional icon  on your desktop, or click Start > All Programs > Agilent > MassHunter Workstation > Mass Profiler Professional > Mass Profiler Professional. b Go to Step 2, Step 3, or Step 4.	MPP begins with the Startup dialog box to help you quickly open a previous project: Open recent project (go to Step 2) Open existing project (go to Step 3) Create new project (go to Step 4) and import an experiment from a previous project
Open recent project: 2 Open a recent project in the Startup dialog box and select the Malaria project from the recent project list.	a Click Open recent project. b Select Malaria from the Select recent project list. c Click OK.	Open recent project opens the project and the experiment(s) that are stored in the project. If your project is not shown in the Select recent project list go to Step 3 and select it from the list of existing projects on your computer. - A project is a container for a collection of experiments. A project can have multiple experiments on different sample types and organisms. - You can now use the operations available in the Workflow Browser.



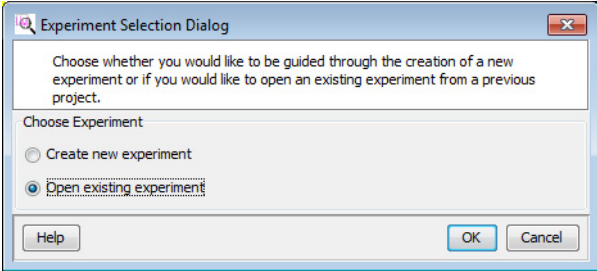
Steps	Detailed Instructions	Comments
Open existing project: 3 Open an existing project in the Startup dialog box and select the Malaria project from the list of all of your existing projects.	a Click Open existing project in the Startup dialog box in Step 2 on page 20 . b Click OK . c Select the Malaria project row from the Open Project dialog box. d Click Open .	• Open existing project allows you to open a project that is no longer displayed in the recent project list. • Open existing project opens the project and the experiment(s) that are stored in the project. • When you have a large number of existing projects on your computer you may have to use the page navigation buttons in the Open Project dialog box. • A selected project is highlighted by a color based on your PC's system settings. • You can now use the operations available in the Workflow Browser.

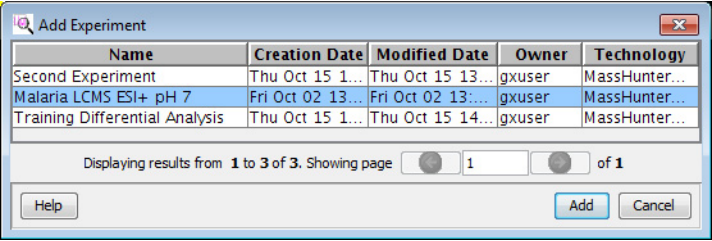


Create new project: 4 Create a new project in the Startup dialog box and import the experiment stored in the Malaria project into the new project.	a Click Create new project in the Startup dialog box in step 2 on page 20 . b Click OK . c Type Malaria Project in Name . d Type Import Experiment in Notes . e Click OK .	• Create new project allows you to import an experiment from an existing project into the new project. • See “Set up a new project” on page 15 for creating a new project and a new experiment.
--	---	---



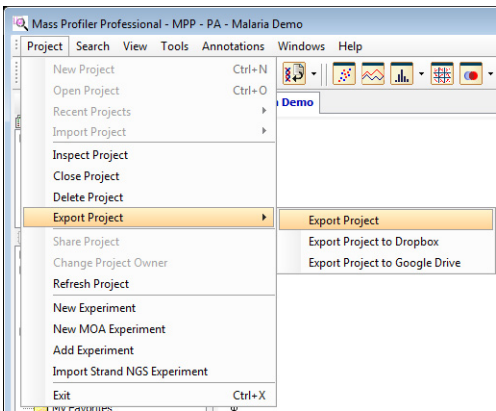
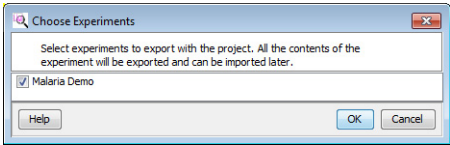
3. Basic Tasks

Steps	Detailed Instructions	Comments
	f Click Open existing experiment in the Experiment Selection Dialog dialog box.	
	 The 'Experiment Selection Dialog' box has a title bar with a magnifying glass icon and a close button. The main text says: 'Choose whether you would like to be guided through the creation of a new experiment or if you would like to open an existing experiment from a previous project.' Below this is a section 'Choose Experiment' with two radio buttons: 'Create new experiment' and 'Open existing experiment'. The 'Open existing experiment' button is selected. At the bottom are 'Help', 'OK', and 'Cancel' buttons.	
	g Select the Malaria LCMS ESI+ pH 7 experiment in the Add Experiment dialog box h Click Add.	• The selected experiment is shown highlighted using a background color. • The <i>Malaria LCMS ESI+ pH 7</i> experiment in the <i>Malaria</i> project is now imported into your new project and appears as if it was opened using Open recent project. • You can now use the operations available in the Workflow Browser.



Save your project

Save your current analysis as a TAR file for archiving, restoration of any future analysis to the current results, sharing the data with a collaborator, or sharing the data with Agilent customer support.

Steps	Detailed Instructions	Comments
1 Export your project to a TAR file.	a Click Project > Export Project.  b Mark the check box next to the experiment you wish to save c Click OK.  d Select or create the file folder. e Type the File name. f Click Save.	

4. Advanced Tasks

The operations available in the Workflow Browser (see [Figure 11](#) on page 13) provide the tools necessary for analyzing features from your mass spectrometry data depending upon the need and aim of the analysis, the experiment design, and the focus of the study. This helps you create different interpretations to carry out the analysis based on the different filtering, normalization, and standard statistical methods.

Follow the procedures in the [Mass Profiler Professional Familiarization Guide](#) (p/n G3835-90026) and [Mass Profiler Professional Application Guide](#) (p/n G3835-90028) (see “Documents” on page 4) for increasingly detailed information to help you use MPP.

The [Agilent Metabolomics Workflow - Discovery Workflow Guide](#) (p/n 5990-7067EN), [Integrated Biology with Agilent Mass Profiler Professional - Workflow Guide Overview](#) (p/n 5991-1910EN), [Class Prediction with Agilent Mass Profiler Professional Workflow Overview](#) (p/n 5991-1912EN), and the [Agilent MassHunter Mass Profiler Professional User Manual](#). (see “Documents” on page 4) provide you with techniques and explanations to perform advanced analysis tasks.

5. Acknowledgment and Citations

BioCyc Pathway/Genome Databases



Includes BioCyc Pathway/Genome databases from the Bioinformatics Research Group at SRI International®, used under license.

<http://www.biocyc.org/>

Citation based on use of BioCyc databases or the Pathway Tools software

If you use BioCyc databases or the Pathway Tools software in your research, cite relevant publications as described on the BioCyc website:

<http://biocyc.org/publications.shtml>

For example, users who publish research results in scientific journals based on use of data from the EcoCyc Pathway/Genome database should cite:

Keseler et al., *Nucleic Acids Research* **39**:D583-90, 2011.

Users who publish research results in scientific journals based on use of data from most other BioCyc Pathway/Genome databases should cite:

Caspi et al., *Nucleic Acids Research* **40**:D742-53, 2012.

KEGG Database



Includes KEGG (Kyoto Encyclopedia of Genes and Genomes) databases developed by Kanehisa Laboratories.

<http://www.genome.jp/kegg/>

Citation based on use of KEGG Database

If you use the KEGG database in your research, cite relevant publications as described on the KEGG website:

<http://www.genome.jp/kegg/kegg1.html>

www.agilent.com

In this book

*The Agilent G3835AA
MassHunter Mass Profiler
Professional Software - Quick
Start Guide* presents the first
steps to use the MassHunter
Mass Profiler Professional
Software.

This guide applies to MassHunter
Mass Profiler Professional
Software 14.0 and higher until
superseded.

© Agilent Technologies, Inc. 2015

Revision A, November 2015



G3835-90025



Agilent Technologies