



Agilent G3835AA MassHunter Mass Profiler Professional Software

Quick Start Guide

Getting Started	4
User Interface	8
Basic Tasks	12
Advanced Tasks	21

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, ICP-MS, and NMR 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.



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 Databases (METLIN, Pesticides, Forensics) 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.
- 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 now can be used with the new, standalone Automated Sample Class Prediction software to enable high-throughput routine classification of unknown samples.

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

MPP Familiarization Guide Agilent G3835AA MassHunter Mass Profiler Professional Software - Familiarization Guide (Agilent publication G3835-90010, Revision A, November 2012)

MPP Application Guide Agilent G3835AA MassHunter Mass Profiler Professional Software - Application Guide (Agilent publication G3835-90011, Revision A, November 2012)

Mass Profiler Professional User Manual Agilent MassHunter Mass Profiler Professional User Manual (Agilent publication, January 2012). You can find a PDF copy of the user manual in the MPP installation folder
C:\Program Files\Agilent\MassHunter\Workstation\Mass Profiler Professional\docs>manual.

Discovery Workflow Guide Agilent Metabolomics Workflow - Discovery Workflow Guide (Agilent publication 5990-7067EN, Revision B, October 2012)

Discovery Workflow Overview Agilent Metabolomics Workflow - Discovery Workflow Overview (Agilent publication 5990-7068EN, Revision B, October 2012)

Qualitative Analysis Familiarization Guide for LC/MS Agilent MassHunter Workstation Software Qualitative Analysis (Agilent publication G3335-90146, Revision A, October 2012)

Qualitative Analysis Familiarization Guide for GC/MS Agilent MassHunter Workstation Software Qualitative Analysis (Agilent publication G3335-90147, Revision A, October 2012)

Quantitative Analysis Familiarization Guide Agilent MassHunter Workstation Software Quantitative Analysis (Agilent publication G3335-90061, Fourth Edition, April 2010)

Training

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


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

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 5).
- 3 Review the following three sections in this *Quick Start Guide*:
“[User Interface](#)” on page 8
“[Basic Tasks](#)” on page 12
“[Advanced Tasks](#)” on page 21
- 4 Review the “[MPP Familiarization Guide](#)” (see “[Documents](#)” on page 3). 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 “[MPP Application Guide](#)” (see “[Documents](#)” on page 3). 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 11).

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, mark your additional products.
- 2 Enter your **Mass Profiler Professional Order ID**. If you have additional software product licenses, enter your additional product Order ID's.
- 3 Click **OK**.

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

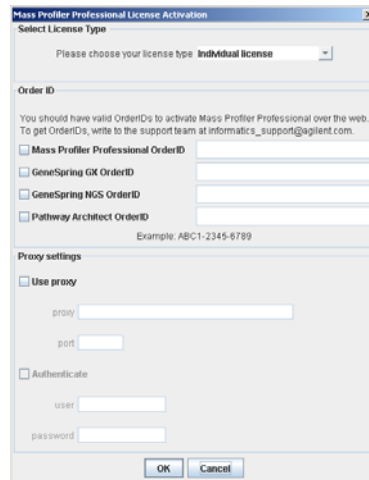


Figure 1 Mass Profiler Professional License Activation dialog box

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 un-install 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.

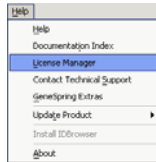


Figure 2 License Manager is available from the Help menu.

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

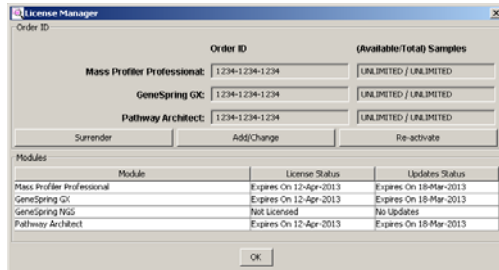


Figure 3 License Manager dialog box

3 Mark **Mass Profiler Professional** and any additional product Order ID's to surrender.

4 Click **OK**.

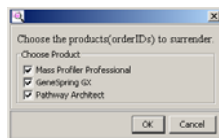


Figure 4 Order ID surrender product selection dialog box

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**.

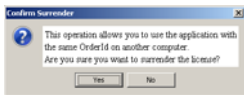


Figure 5 Confirm Surrender dialog box

- 6 Click **OK** in the **Automatic Surrender Operation Successful** dialog box. Because your license is no longer active on the current computer, MPP software immediately closes.

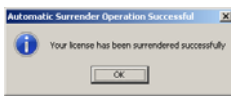


Figure 6 Automatic Surrender Operation Successful dialog box

- 7 Activate your license Order ID on the new computer (see "[Activating a License](#)" on page 5).
- 8 *(Optional)* Click **Start > All Programs > Agilent > MassHunter Workstation > Mass Profiler Professional > Uninstall** to remove MPP from the current computer.

User Interface

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*.

Main Functional Areas

The main functional areas of Mass Profiler Professional are shown in [Figure 9](#) on page 9.

The main MPP window consists of four parts: (1) the Menu Bar, (2) the Toolbar, (3) the Display Pane, and (4) the Status Bar.

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

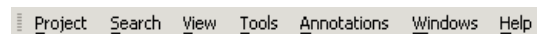


Figure 7 Menu bar

2. Toolbar The toolbar ([Figure 8](#)) 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 and Add experiment

Entity List section: Create entity list from selection, Inspect selected entity, and Import entity list from file

Plots, Statistics, and Pathways section: Scatter plot, Profile plot, Histogram plot, Matrix plot, Venn diagram plot, Box-and-whisker plot, Heat map plot, Data spreadsheet, Summary statistics, Create new pathway, and Select data source for plots.

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



Figure 8 Toolbar

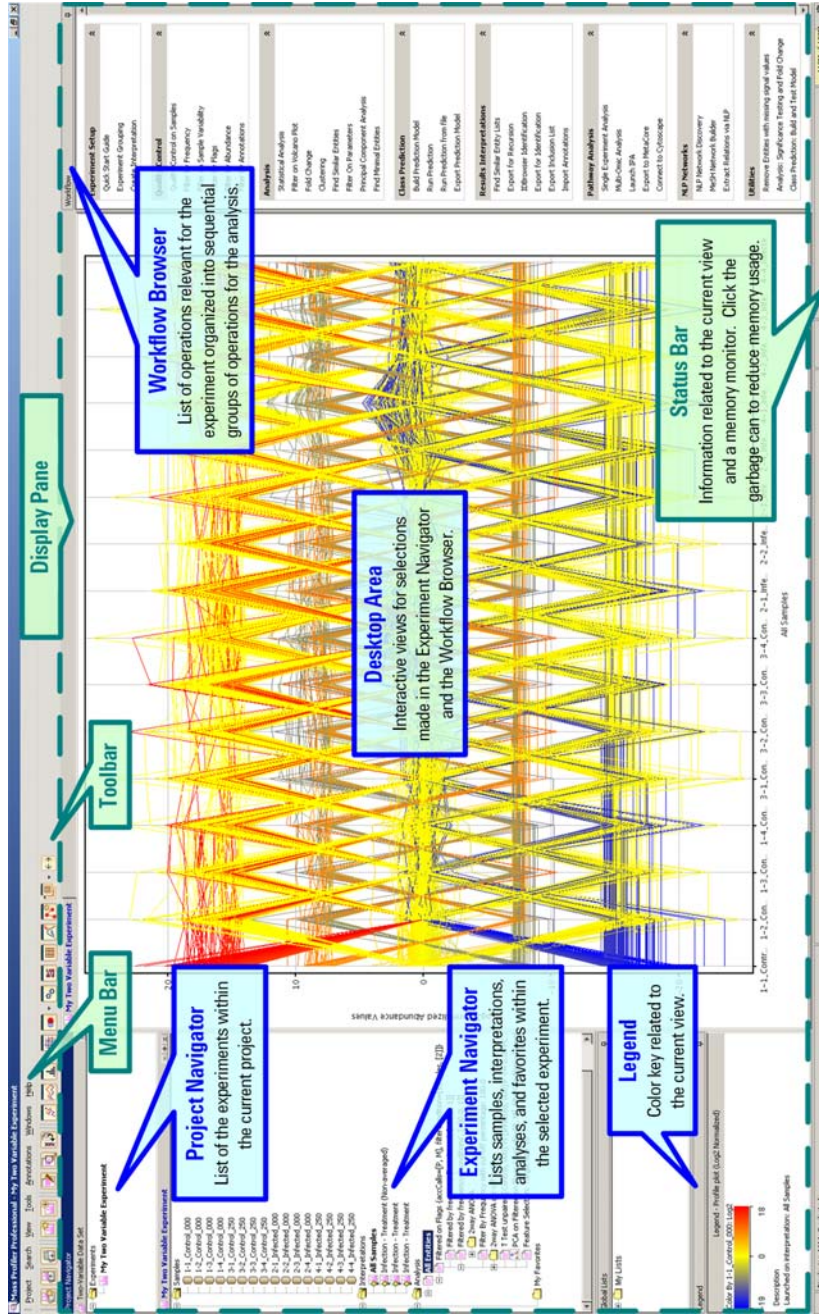


Figure 9 The main functional areas of the MPP software

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

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

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](#)) 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.

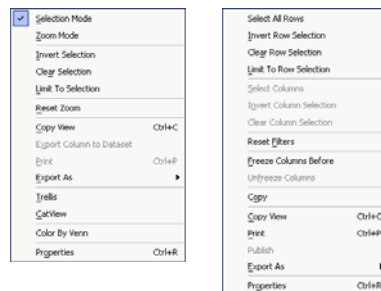


Figure 10 The right-click available options within the desktop view automatically adjust to the type of view.

Workflow Browser

The Workflow Browser (Figure 11, left) 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.

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 described in the Desktop area.

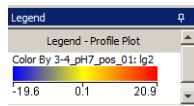


Figure 12 Legend

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.

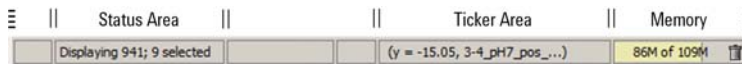


Figure 13 Status Bar

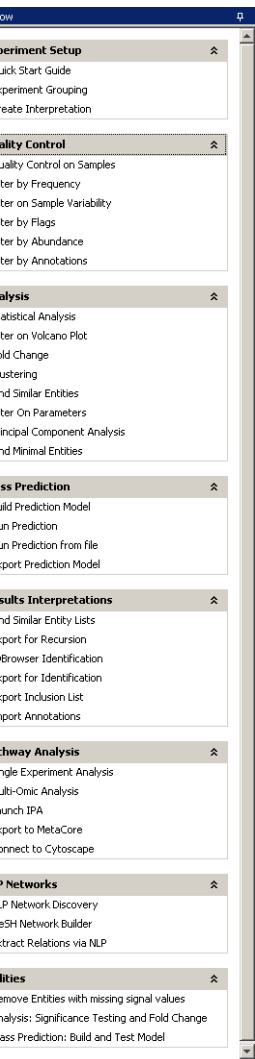


Figure 11 Workflow Browser operations



Basic Tasks

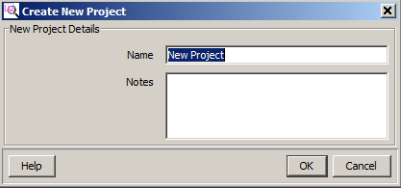
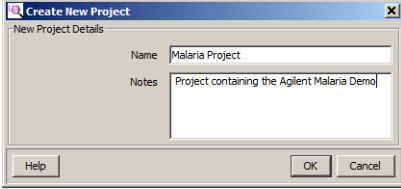


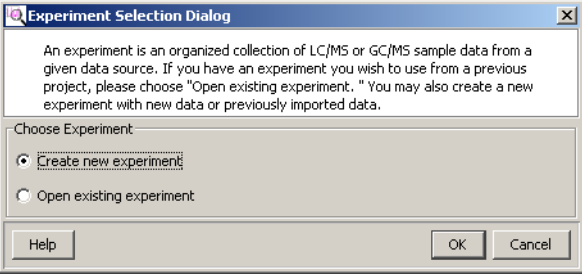
Set up a new project

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.	<p>a Click the Mass Profiler Professional icon  on your desktop, or click Start > All Programs > Agilent > MassHunter Workstation > Mass Profiler Professional > Mass Profiler Professional.</p>	<ul style="list-style-type: none">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 17).
2 Create a new project from the Startup dialog box.	<p>a Click Create new project. b Click OK.</p>	<ul style="list-style-type: none">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.

Steps	Detailed Instructions	Comments
<p>3 Type descriptive information in the New Project Details section of the Create New Project dialog box.</p>	<p>a Type <code>Malaria Project</code> in Name. b Type <code>Project containing the Agilent Malaria Demo</code> in Notes. c Click OK.</p>	<ul style="list-style-type: none"> 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.
<div style="display: flex; align-items: center; justify-content: center;">  >  </div>		
<p>4 Create a new experiment in the Experiment Selection Dialog dialog box.</p>	<p>a Click Create new experiment. b Click OK.</p>	<ul style="list-style-type: none"> 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 17.
<div style="display: flex; align-items: center; justify-content: center;">  </div>		

Steps	Detailed Instructions	Comments
5 Type and select information that guides the experiment creation in the New Experiment dialog box.	<p>a Type the descriptive name <code>Malaria demo</code> in Experiment name. This name may be different from the project name.</p> <p>b Select Mass Profiler Professional for Analysis type.</p> <p>c Select Unidentified for the Experiment type.</p> <p>d Select Analysis: Significance Testing and Fold Change for Workflow type.</p> <p>e Type <code>Malaria LCMS ESI+ pH 7</code> in Experiment notes.</p> <p>f Click OK.</p>	<ul style="list-style-type: none"> • Experiment type determines how Mass Profiler Professional manages the data. • Unidentified is the proper Experiment type selection when the compounds have only been identified by their molecular features of neutral mass and retention time. • Identified is the proper Experiment type selection when the compounds have been identified by compound, formula, and/or CAS number. • 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 MassHunter Qualitative Analysis has been previously used to identify some of the compound features.



- 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 “[MPP Familiarization Guide](#)” (see “[Documents](#)” on page 3) guides you through importing and organizing the *Malaria Demo* data.

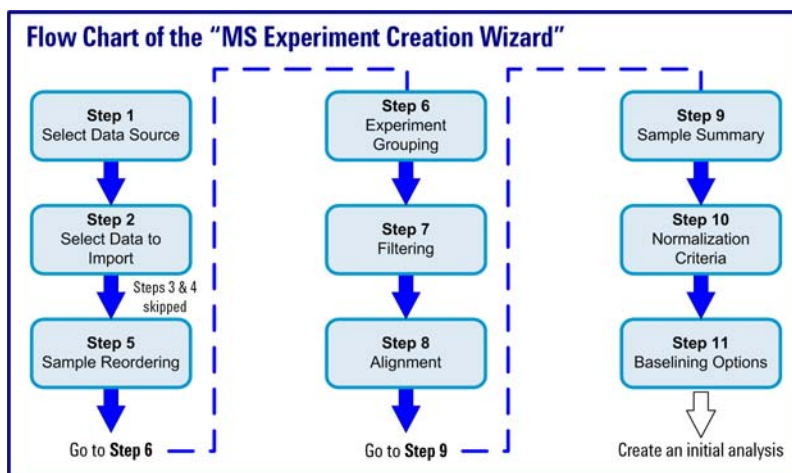


Figure 14 “MS Experiment Creation Wizard”

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. The “MPP Familiarization Guide” (see “Documents” on page 3) guides you through an initial differential analysis using the *Malaria Demo* data.

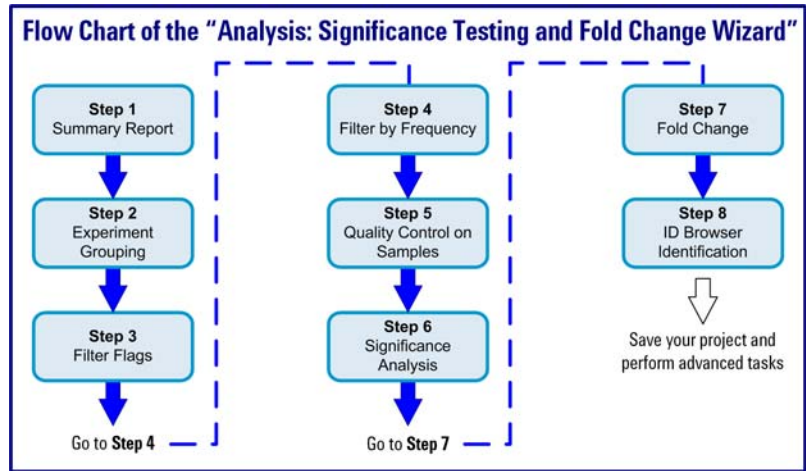



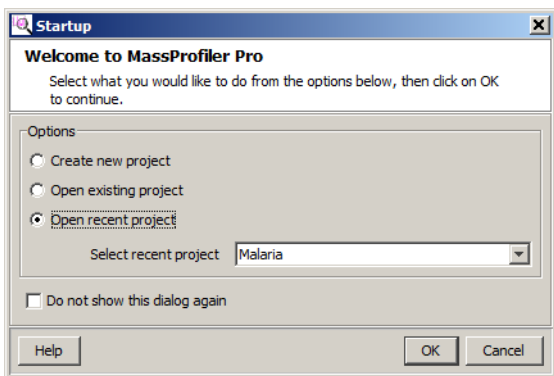
Figure 15 “Analysis: Significance Testing and Fold Change Wizard”

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 12 and then refer to the “[MPP Familiarization Guide](#)” (see “[Documents](#)” on page 3).

Steps	Detailed Instructions	Comments
1 Start Mass Profiler Professional.	<p>a Click the Mass Profiler Professional icon  on your desktop, or click Start > All Programs > Agilent > MassHunter Workstation > Mass Profiler Professional > Mass Profiler Professional.</p> <p>b Go to Step 2, Step 3, or Step 4.</p>	<ul style="list-style-type: none"> 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
<p>Open recent project:</p> <p>2 Open a recent project in the Startup dialog box and select the Malaria project from the recent project list.</p>	<p>a Click Open recent project.</p> <p>b Select Malaria from the Select recent project list.</p> <p>c Click OK.</p>	<ul style="list-style-type: none"> 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 may now use the operations available in the Workflow Browser.



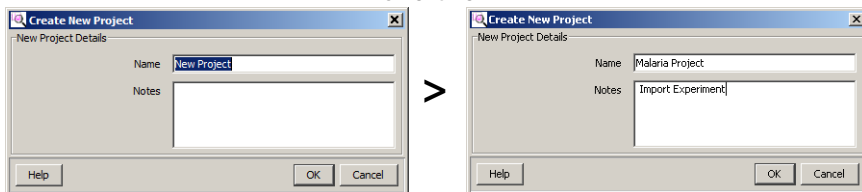
Steps	Detailed Instructions	Comments
-------	-----------------------	----------

Create new project:

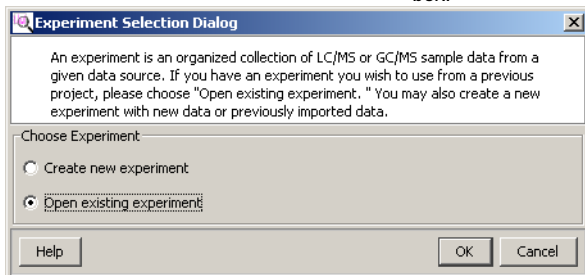
4 Create a new project in the **Startup** dialog box and import the experiment stored in the Malaria project into the new project.

- a Click **Create new project** in the **Startup** dialog box in step 2 on page 17.
- 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 12 for creating a new project and a new experiment.

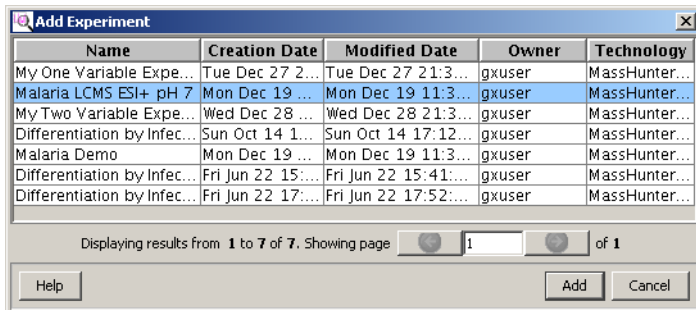


f Click **Open existing experiment** in the **Experiment Selection Dialog** dialog box.



- 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 *Malaria LCMS ESI+ pH 7* experiment in the *Malaria* project is now imported into your new project and appears as if it was opened using **Open recent project**.
- You may now use the operations available in the Workflow Browser.

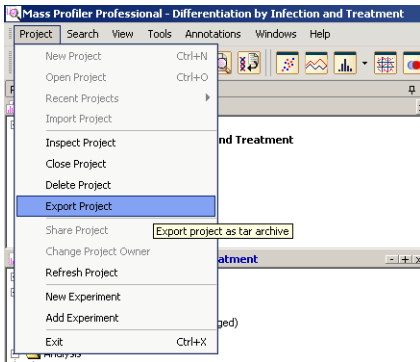


Save your project

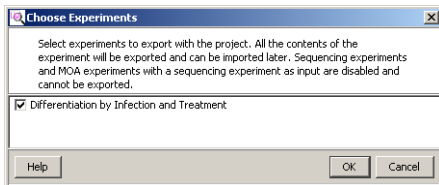
Save your current analysis as a TAR file for archiving, restoration of any future analysis to the current results, sharing the data with a collaborator, or sharing the data with Agilent customer support.

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**.

Advanced Tasks

The operations available in the Workflow Browser (see [Figure 11](#) on page 11) 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.

It is recommended that you follow the procedures in the “[MPP Familiarization Guide](#)” and “[MPP Application Guide](#)” (see “[Documents](#)” on page 3) for increasingly detailed information to help you use MPP.

The “[Discovery Workflow Guide](#)” and the “[Mass Profiler Professional User Manual](#)” (see “[Documents](#)” on page 3) provide you with techniques and explanations to perform advanced analysis tasks.

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

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.

In some cases, BioCyc Pathway/Genome databases are described by other specific publications that can be found by selecting the database and then going to the Summary Statistics pages under the Tools menu. The resulting page sometimes contains a citation for that database.

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.

© Agilent Technologies, Inc. 2012

Revision A, November 2012



G3835-90009



Agilent Technologies