



Agilent MassHunter GC/MS System

Quick Start Guide



Notices

© 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.

Manual Part Number

G7006-90014

Edition

First Edition, October 2024

Printed in USA

Agilent Technologies, Inc.
5301 Stevens Creek Boulevard
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

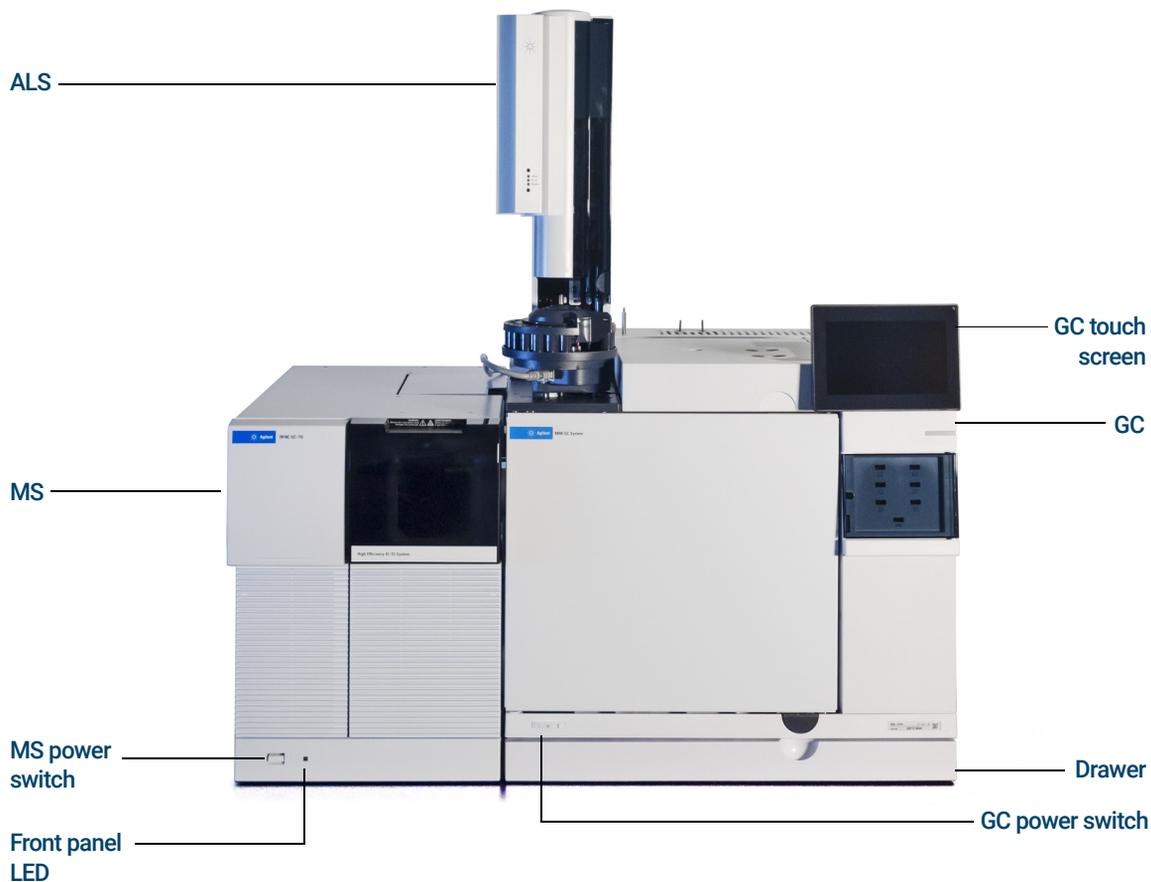
A CAUTION notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a CAUTION notice until the indicated conditions are fully understood 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.

The Agilent GC/MS System

Below is a typical Gas Chromatograph and Mass Spectrometer (GC/MS) hardware configuration.



The Agilent MassHunter Software Suite

The MassHunter software suite offers dedicated software modules for compound quantitation and confirmation, target compound screening, degradant and unknowns identification, characterization of biomolecules, and protein and metabolite identification.

To open a module, select the corresponding desktop icon:



Data Acquisition



Report Builder



Quantitative Analysis



Library Editor



Qualitative Analysis



Unknowns Analysis



OpenLab Control Panel



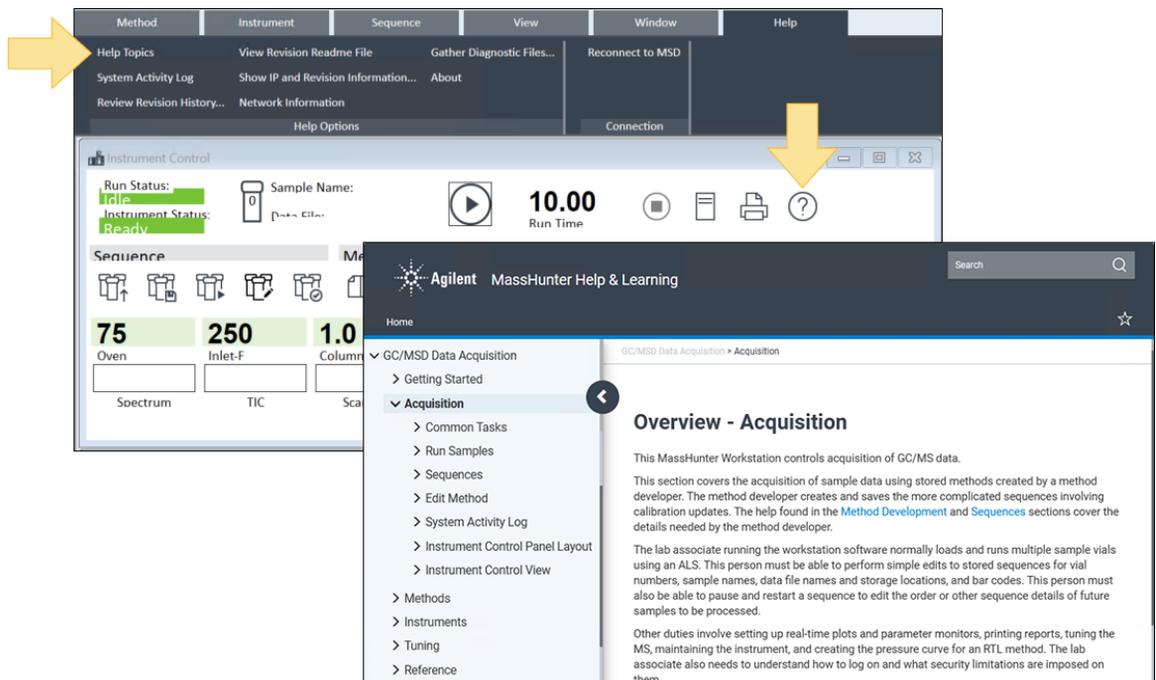
Parts Finder

MassHunter Online Help

An extensive library of software information and tutorials about instrument control, data acquisition, data analysis, methods, sequencing, tuning, troubleshooting, and how to use system commands and variables is available through the online Help files.

To access the online Help from your MassHunter software, do one of the following:

- Press F1.
- Select the  Help icon.
- On the **Help** tab, select **Help Topics**.





Data Acquisition Software

From MassHunter Data Acquisition, you can:

- Start and stop the instrument.
- Download instrument settings in real time.
- Load, edit, create, or run a method.
- Load, edit, create, check, or run a sequence.
- Evaluate the MS parameters.
- Autotune the MS parameters.
- Monitor the real-time conditions of the instrument parameters.
- View the instrument logs.
- View and print the real-time plot for chromatograms and instrument spectra.
- Select and label the total ion chromatograms (TIC) or extracted ion chromatograms (EIC) that you want to appear in the real-time plot.

Select a tab to view additional windows and options.



Hover over any icon or button to display a tooltip identifying what each tool can do.



| Method | Instrument | Sequence | View | Window | Help |
|-------------------|-------------------------|---|------|----------------------------------|-----------------|
| Load Method... | Run Method... | Set/Change Standby Method <None>... | | Edit Entire Method... | 1 SampleA-RTL.m |
| Save Method | Print Method... | Import Method Components From Another Method... | | Additional Method Information... | 2 |
| Save Method As... | View Method Audit Trail | | | Edit Method | 3 |

Method Actions

Instrument Control

Run Status: 10.00 Run Time

Sample Name: [...]

Instrument Status: [...]

Sequence

Oven: 75

Inlet-F: 250

Column-1 Flow: 1.0

Guard Chio: 0

MS Transfer Line: 250

MS Source (°C): 230

MS Quad (°C): 150

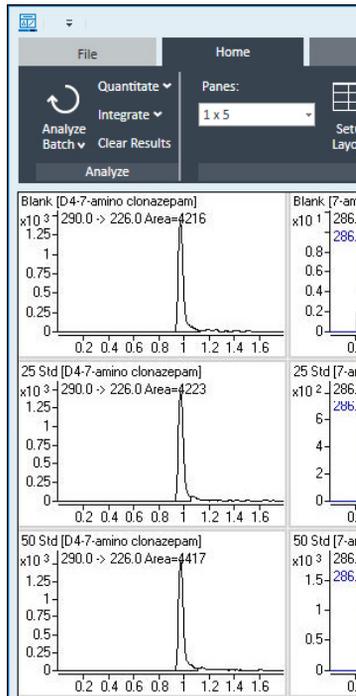
Spectrum

TIC

Scan 1-1



Quantitative Analysis Software



The MassHunter Quantitative Analysis software will quantify very low amounts of sample.

From one screen in MassHunter Quantitative Analysis, you can:

- Import information directly from the acquisition method.
- Use the curve-fit assistant to test all fits and statistics on the calibration curve quality.
- Use the Compounds at a Glance feature to review all the chromatograms in your batch at once, side-by-side.
- Quickly scan the batch table and see your quality metrics such as high outliers (in red) or low outliers (in blue).

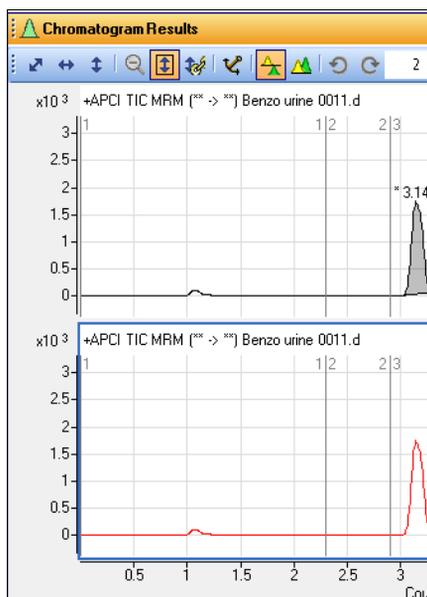
You may access familiarization guides for this software through the online Help.

Press F1 to access context sensitive online Help.

| Batch Table | | | | | | | | | | |
|---|----------------------------|---------------------------------------|--------------------------|-------------|-------------|----------|--------------|--------------------------|-----------------|-------|
| Sample: 7-amino clonazepam | | Sample Type: <All> | | | | | | | | |
| 7-amino c... | 7-amino clonazepam Results | | | | | | Qualifier... | | D 4-7-amino ... | |
| Exp. Conc. | RT | Resp. | MI | Calc. Conc. | Final Conc. | Accuracy | Ratio | MI | RT | Resp. |
| 25.0000 | 0.982 | 2400 | <input type="checkbox"/> | 29.0230 | 29.0230 | 116.1 | 79.1 | <input type="checkbox"/> | 0.974 | 4223 |
| 50.0000 | 0.982 | 5192 | <input type="checkbox"/> | 48.6966 | 48.6966 | 97.4 | 79.2 | <input type="checkbox"/> | 0.974 | 4417 |
| 125.0000 | 0.982 | 12941 | <input type="checkbox"/> | 111.0894 | 111.0894 | 88.9 | 85.4 | <input type="checkbox"/> | 0.973 | 4173 |
| 250.0000 | 0.984 | 28338 | <input type="checkbox"/> | 227.0264 | 227.0264 | 90.8 | 82.6 | <input type="checkbox"/> | 0.977 | 4243 |
| 500.0000 | 0.983 | 60068 | <input type="checkbox"/> | 534.1646 | 534.1646 | 106.8 | 74.5 | <input type="checkbox"/> | 0.975 | 3717 |
| | | | <input type="checkbox"/> | | | | | <input type="checkbox"/> | 0.972 | 4216 |
| | | | <input type="checkbox"/> | | | | | <input type="checkbox"/> | 0.976 | 4599 |
| 189.0000 | 0.982 | 17646 | <input type="checkbox"/> | 145.8304 | 145.8304 | 77.2 | 75.8 | <input type="checkbox"/> | 0.974 | 4228 |



Qualitative Analysis Software



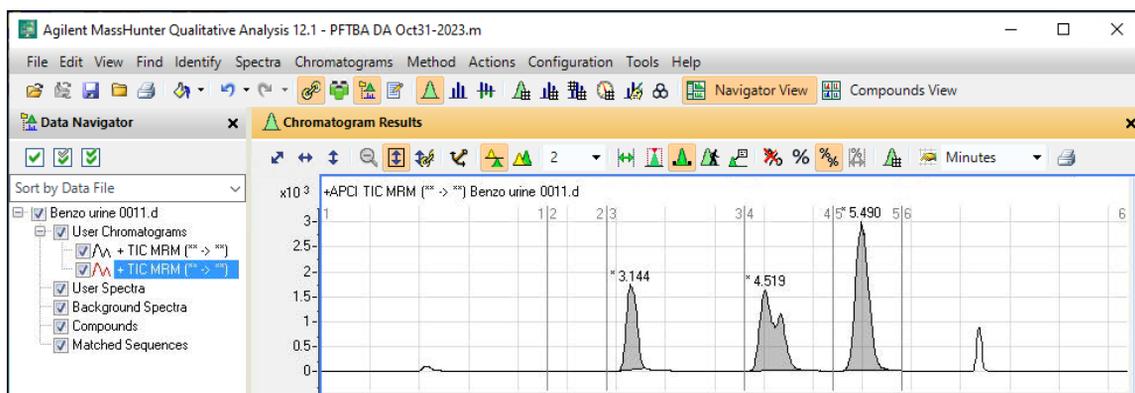
The MassHunter Qualitative Analysis software uses advanced data mining and processing tools, including integrated deconvolution algorithms, to rapidly and accurately find all detectable compounds in your samples and easily confirm targets or identify unknowns.

This software presents large amounts of data for review in one central location. From one screen, you can:

- Extract chromatograms
- View and extract peak spectra
- Subtract background
- Integrate the chromatogram
- Find compounds and generate formulas

You can also set up methods to automatically do the tasks in the above list, as well as others, when you open the data files.

You may access familiarization guides for this software through the online Help. Press F1 to access context sensitive online Help.





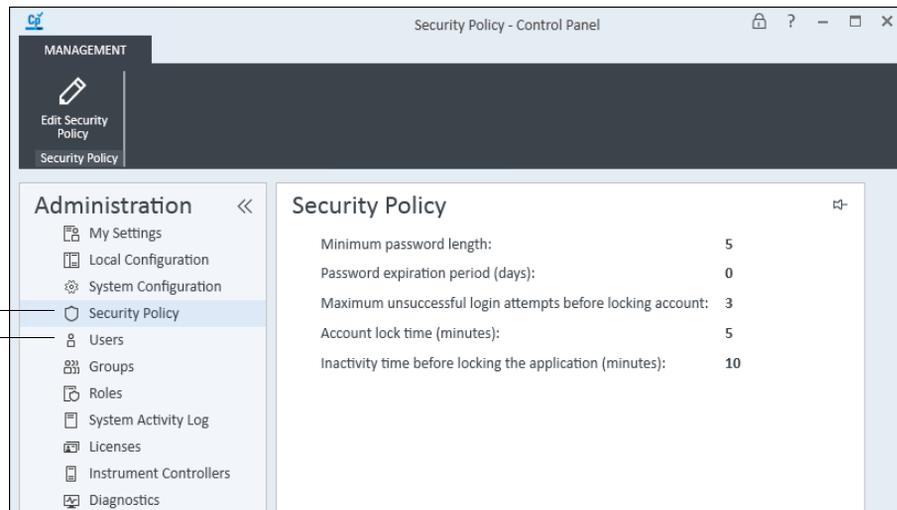
OpenLab Control Panel

The Control Panel is the administrative and management center of MassHunter, offering:

- Full instrument status information for your entire laboratory
- Central configuration and administration of users, instruments, and security settings
- Full system documentation and built-in reports.

Select **Security Policy** to set the login Security Policy and control your users' access to the system.

Select **Users** to add users to your system and set their privileges.



For details on the Control Panel features, press F1 to see the online Help.



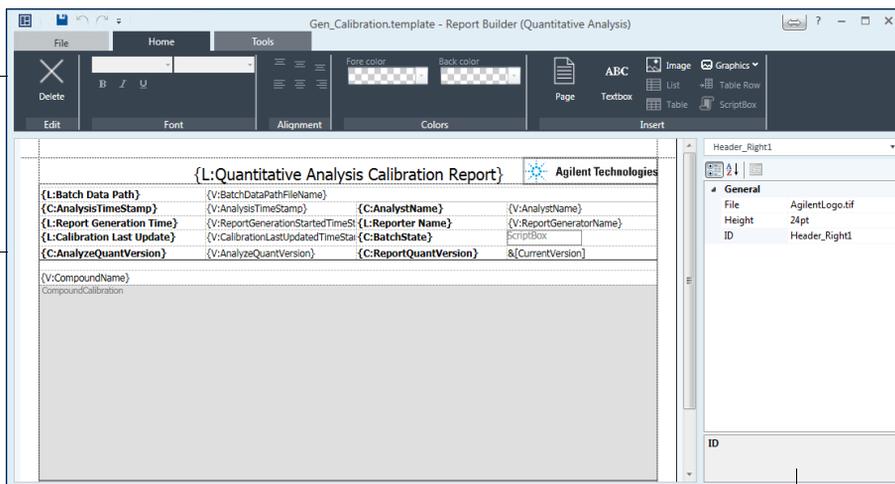
Report Builder

The Report Builder software is installed with MassHunter Quantitative Analysis. The software provides PDF templates for MassHunter Quantitative Analysis and MassHunter Unknowns Analysis. Each of these templates can be customized for your specific application in the Report Builder.

You may access familiarization guides for this software through online Help.

Use the toolbar on the **Home** tab to edit fonts, colors, and alignment and insert various elements into the template.

Use the editing window to preview the **report template**.



Use the **properties** pane to edit details for each element.



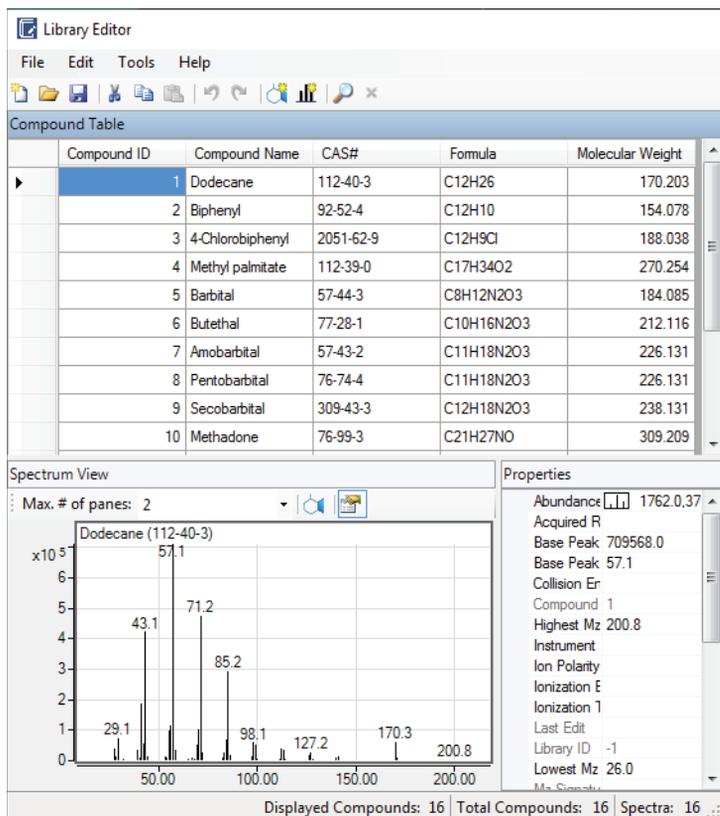
Library Editor Software

The MassHunter Library Editor software customizes libraries used to identify compounds and components in the MassHunter Qualitative and Quantitative Data Analysis software.

From one screen in MassHunter Library Editor, you can:

- Add compounds to a library
- Copy spectral data from your data analysis software to a library
- Add a structure to your custom library
- Import JCAMP files into a custom library
- Create a reference pattern library

You may access familiarization guides for this software through the online Help.





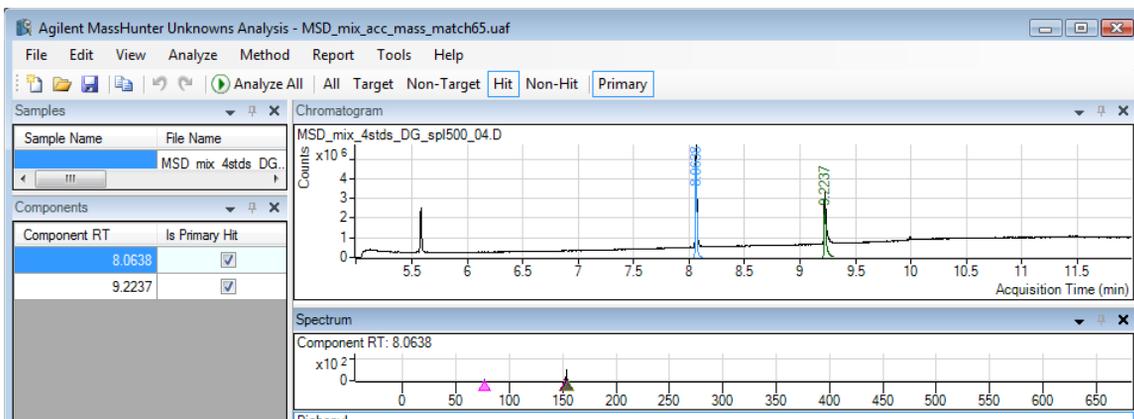
Unknowns Analysis Software

The MassHunter Unknowns Analysis software identifies compounds that may be present in a sample batch beyond those that were found in MassHunter Quantitative Analysis.

From one screen in MassHunter Unknowns Analysis, you can:

- Import data directly from MassHunter Quantitative Analysis.
- Use optimized deconvolution and library match algorithms to identify target and non-target hits.
- View chromatograms, component lists, spectrums, ion peaks, and molecular structures side-by-side.
- Easily review library match and peak shape metrics of unknowns components in table and spectrum formats.

You may access familiarization guides for this software through the online Help.





Parts Finder Tool

Included with your system is the Agilent Parts Finder tool. This interactive software features an intuitive image-based search capability for quick identification of needed parts and supplies for your Agilent instruments.

Using Parts Finder, rather than thumbing through a paper catalog, you can select nested photos to quickly locate the parts that you need.



Parts Finder is installed on your PC along with MassHunter Workstation.

MS Hardware Library

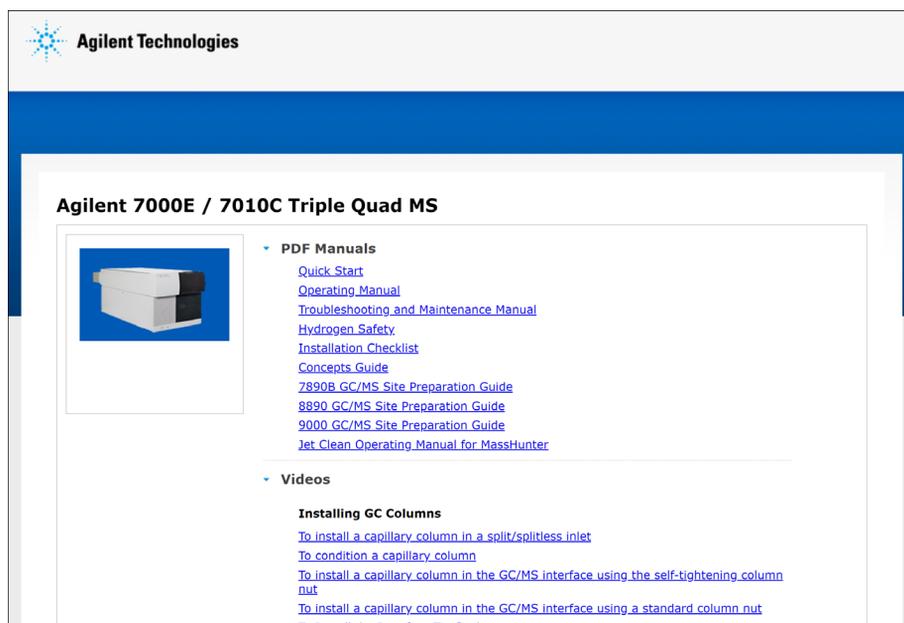
Use the Agilent GC/MS Hardware User Information USB to install the full library of Learning Products with manuals (in PDF format) and videos (in MP4 format) that explain how to install, operate, maintain, and troubleshoot your system.



To install your MS Hardware Library, insert the memory stick into a USB port and follow the prompts. This can be installed by anyone who has authority to copy information onto the receiving computer.



Once installed, an icon appears on your desktop for quick access to your material. To open a hardware library, select the desktop icon.



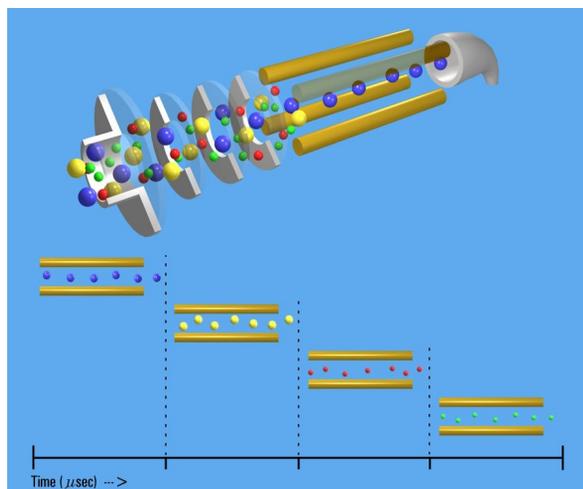
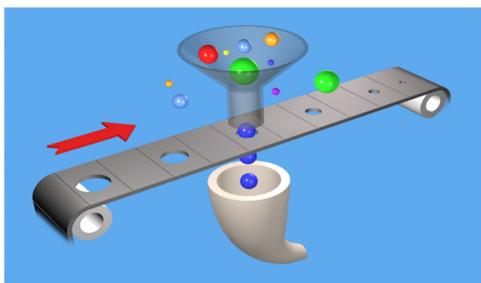
The screenshot displays the Agilent Technologies website interface. At the top left, the Agilent Technologies logo and name are visible. Below this, a blue header bar spans the width of the page. The main content area features a white box with a blue border. On the left side of this box is a photograph of the Agilent 7000E / 7010C Triple Quad MS instrument. To the right of the image, the text "Agilent 7000E / 7010C Triple Quad MS" is displayed. Below the image and title, there are two sections: "PDF Manuals" and "Videos". The "PDF Manuals" section lists several links: Quick Start, Operating Manual, Troubleshooting and Maintenance Manual, Hydrogen Safety, Installation Checklist, Concepts Guide, 7890B GC/MS Site Preparation Guide, 8890 GC/MS Site Preparation Guide, 9000 GC/MS Site Preparation Guide, and Jet Clean Operating Manual for MassHunter. The "Videos" section is titled "Installing GC Columns" and lists links for: To install a capillary column in a split/splitless inlet, To condition a capillary column, To install a capillary column in the GC/MS interface using the self-tightening column nut, To install a capillary column in the GC/MS interface using a standard column nut, and To Install the Interface Tin Seal.

Concepts Guide

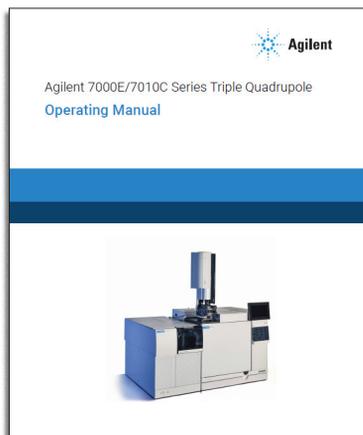


To learn how your Agilent GC/MS system works in conjunction with the MassHunter software, take a look at the concepts guide included with your Hardware Library. This document includes:

- Conceptual models
- Illustrations
- Screen captures
- An overview of how the mass spec works
- Concepts behind the design of MassHunter Data Acquisition
- The theory behind chemical ionization in GC/MS



Operating Manual

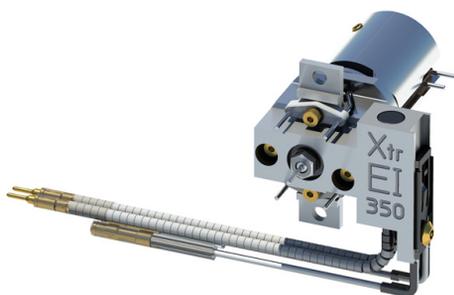


Your Agilent GC/MS system is extremely powerful and flexible. Refer to the operating manual included with your Hardware Library for extensive information on operating the instrument and performing routine maintenance on it. Included are step-by-step instructions and videos on installing GC columns, operating in electron impact (EI) mode, general maintenance procedures, and more.



Any topic with the camera icon is linked to a video. Watch these videos to learn

more.



Troubleshooting and Maintenance Manual

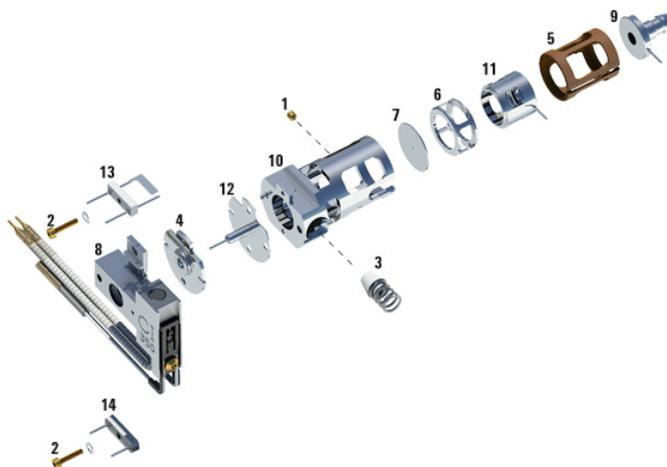


The troubleshooting and maintenance manual included in your Hardware Library includes step-by-step instructions along with videos, IPBs, and annotated high-resolution photos to help you identify the cause of problems in your system. Detailed procedures are also included for completing common maintenance tasks, sections on the components of the GC/MS vacuum system and analyzer system, and a listing of parts.

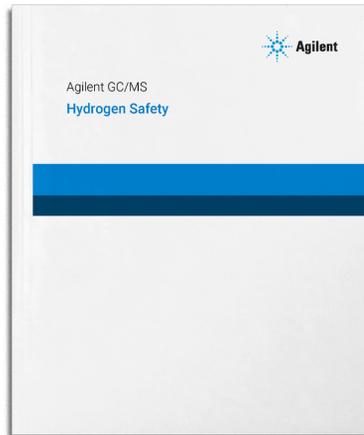


Any topic with the camera icon is linked to a video. Watch these videos to learn

more.



Hydrogen Safety Manual

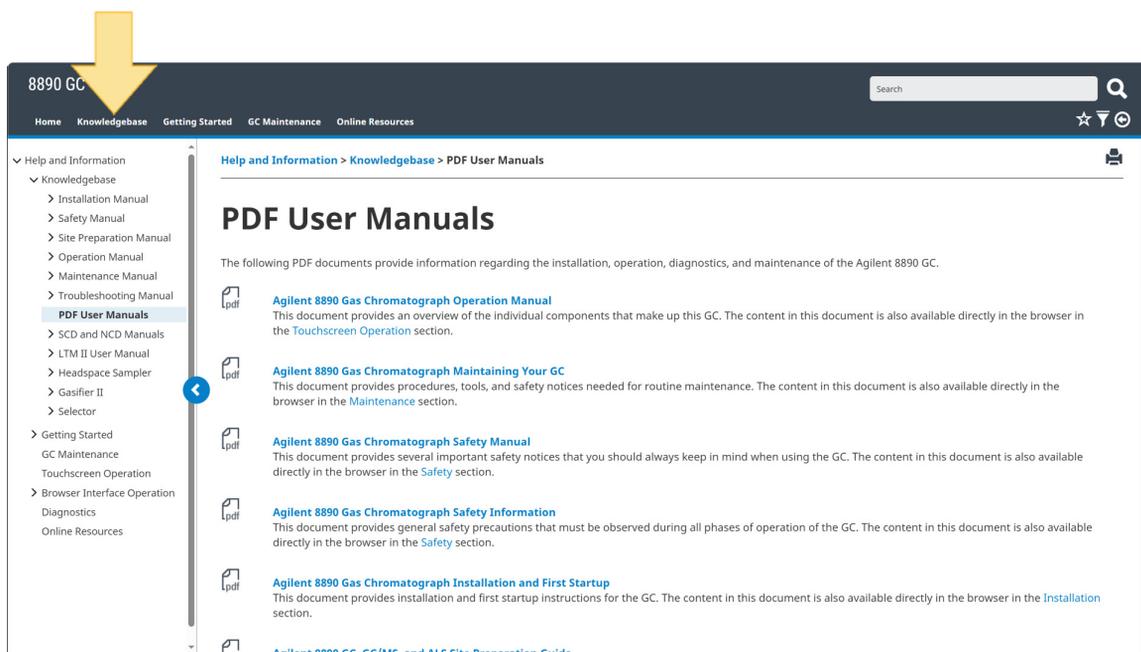


Hydrogen presents a number of dangers. Some are general, others are unique to GC or GC/MS operation. Review the hydrogen safety manual included in your Hardware Library for details on the use of hydrogen with the GC/MS.

GC Hardware Library

Additional information for your GC is also available.

- 1 Open a browser on a computer or other device that shares the same gateway as the GC.
- 2 Enter this URL: `http://xxx.xxx.xxx.xxx/info`, where `xxx.xxx.xxx.xxx` is the IP address or host name of the GC. For example, enter `http://10.1.1.101/info`.
- 3 When the GC Help & Information home screen opens, select **Knowledgebase**.



The screenshot displays the Agilent 8890 GC Knowledgebase interface. A yellow arrow points to the 'Knowledgebase' tab in the top navigation bar. The left sidebar shows a tree view with 'PDF User Manuals' selected. The main content area is titled 'PDF User Manuals' and lists five PDF documents with brief descriptions:

- Agilent 8890 Gas Chromatograph Operation Manual**: This document provides an overview of the individual components that make up this GC. The content in this document is also available directly in the browser in the [Touchscreen Operation](#) section.
- Agilent 8890 Gas Chromatograph Maintaining Your GC**: This document provides procedures, tools, and safety notices needed for routine maintenance. The content in this document is also available directly in the browser in the [Maintenance](#) section.
- Agilent 8890 Gas Chromatograph Safety Manual**: This document provides several important safety notices that you should always keep in mind when using the GC. The content in this document is also available directly in the browser in the [Safety](#) section.
- Agilent 8890 Gas Chromatograph Safety Information**: This document provides general safety precautions that must be observed during all phases of operation of the GC. The content in this document is also available directly in the browser in the [Safety](#) section.
- Agilent 8890 Gas Chromatograph Installation and First Startup**: This document provides installation and first startup instructions for the GC. The content in this document is also available directly in the browser in the [Installation](#) section.
- Agilent 8890 GC GC/MS and ALS Site Preparation Guide**

www.agilent.com

© Agilent Technologies, Inc. 2024

First Edition, October 2024



G7006-90014

