

MS Workstation 7.0.1 (2/29/2012)

Contents

1. New Features.....	1
1.1 Hardware.....	1
1.2 Software.....	1
2. Software Support.....	1
3. Computer Requirements.....	2
4. Modules Supported.....	2
5. Modules Not Supported.....	2
6. Software Supported.....	2
7. Software Not Supported.....	3
8. Known Issues.....	3
9. Injecting Manually, with a CombiPal, or other Third-Party Device.....	3
10. CombiPal SPME Fiber Exposure.....	3

1. New Features

This version of MS Workstation provides support for the MantaRay version of the 7890A Driver on MS Workstation, in addition to, defect fixes.

1.1 Hardware

No new hardware features were added for this release.

1.2 Software

The Dual/Multi Injection and Overlapped Samples features are not supported in MS Workstation 7.0.1.

In addition, due to a 19 character limited for the Sample Name Field in the Sample List, any Barcode Symbologies that have more than 19 characters are not supported (Code 128 (20 characters), Code 128 (40 characters), and Interleaved 2 of 5 (31 characters)).

2. Software Support

For Software Support, please use the following contact information:

US & Canada: 800-227-9770

Worldwide: <http://www.chem.agilent.com/en-us/ContactUS/Pages/ContactUs.aspx>

3. Computer Requirements

The following minimum computer requirements are recommended:

- 1) Intel® Core™2 Duo Processor (or higher)
- 2) Video supporting 1280 x 1024 x 256 resolution or greater
- 3) 32-bit color is recommended
- 4) 2 GB RAM (or more)
- 5) 100 GB Disk (or larger)
- 6) CD/DVD-ROM drive
- 7) Windows 7 Professional or XP

4. Modules Supported

The following modules are supported:

- MS: 2000 MS, 200-MS (210/220/225), 240-MS/4000 MS, 500-MS, 310/320/525 LC/MS
- GC: 450-GC, 431-GC, 3800 GC, 3900 GC, CombiPal
- LC: 212-LC, 460-LC, 8X0 MIB, ProStar 220/230/240, 325, 335, 363, 400, 410, 420, 430, 500, ProStar/Dynamax, Mistral Column Oven (PS 510), CombiPal

5. Modules Not Supported

Support for the following modules was eliminated in this release (or previously):

- MS: 1200 GC/MS, 1200L LC/MS, 300 GC/MS, 920 FTMS
- GC: 39XL GC, 3400/3600 GC, 8200 AS
- LC: 216-LC, ProStar 310, 320, 520, 9001/9002/9010/9012, 9050, 9100, 9200, 9300, ADC Board, AI-200.

6. Software Supported

- Windows XP and 7 Professional (32-bit only)
- EnviroPro, MultiCompound, ToxProPlus, and Dioxins Reporting
- G1033A NIST 11 MS Library
- G1035C Wiley 9th Edition/NIST 11 MS Library
- G1033A Nist 08 MS Library
- G1035B Wiley 9th/NIST 08 Mass Spectral Database
- NIST 08 Mass Spectral Database (Varian P/N: 391082491)
- Wiley 9th Mass Spectral Database (Varian P/N: 391053501 or 391093501)
- Other NIST-formatted libraries
- Access Control and Audit Trail (21CFR11 Support)

7. Software Not Supported

- Any other version of Windows (e.g., Vista, Win 7 64-bit, 2000, NT, ME, 98, 95)
- MS to SMS and DAT to XMS Data File Conversion programs
- QuickAccess Application
- StarFinder, StarWriter, PolyView, and Aurora applications
- LBR2NIST and PalDriv utilities

A Demo version of NIST 08 Library and MS Search Program will be installed as part of the initial MS Workstation installation if the NIST Library is not already installed.

8. Known Issues

Click here ([Software Status Bulletin](#)) for the list of corrected and outstanding issues as of the release date of this version of the software.

9. Injecting Manually, with a CombiPal, or other Third-Party Device

Operation of the CombiPal with a 7890A GC is supported in the MS Workstation 7.0.1 release. Manual, AutoMix, and Automatic Injection Modes (Sample List Injection Mode Options) are supported in the current release. The “No Injection” Injection Mode is NOT supported.

For CombiPal instruments that contain a “Prepare” Object as one of the SyncOut Signals, the GC will be sent a Pre-Run signal just prior to the injection of Sample. For other AutoSamplers, there is a Pre-Run button in the 7890A Module Window in System Control, which allows users to set a specific delay between the start of the AutoSampler Prep actions and entry to Pre-Run on the GC. This will delay the GC from going into the Pre-Run State for the specified amount of time and save on gas usage for long incubation periods or similar prep actions.

Manual Injections are also supported in cases where no AutoSampler is attached to the 7890A GC. If the 7890A is NOT in the AutoStart box on the System Control Configuration Screen, the Pre-Run button can be used to set a delay. The GC will be triggered to enter the Pre-Run State at the end of the specified delay time. If the 7890A is in the AutoStart box on the System Control Configuration Screen, the 7890A will be automatically triggered to go into the Pre-Run State when preparing to do a run. Any delay that is set through the pre-run button on the 7890A Module Window in System Control will be ignored. The user will need to press the Start button on the Front Panel of the 7890A GC (after the Pre-Run is completed) as the user is making the manual injection.

In cases where the 7693 is attached to the 7890A, the Pre-Run button is not enabled (or used) in the 7890A Module Window of System Control.

10. CombiPal SPME Fiber Exposure

In order to adjust the length of the SPME fiber exposed, an entry needs to be added to the “System Control” section of the `\AgilentWS\System\STAR32.ini` file as shown:

[System Control]

CPAL_INJ_FIBER_EXPOSE=XX

The degree of exposure can be adjusted by varying the XX value between 10 and 30. If the entry is missing, the default value is 15.

To have a record of the Vial Penetration and Fiber Exposure values in the Message Log, the following line should also be added to the System Control section:

CPAL_FIBER=1