



Agilent Headspace Driver Release Notes

[B.01.05 Release Note](#)

November 22, 2015

Agilent Headspace Firmware Interoperability and Support Statements

Agilent releases GC firmware updates independently of software releases.

Agilent Headspace instrument driver revisions have been designed to be backward compatible to the installed instrument base.

Agilent recommends always using the latest module firmware revision in order to provide the highest level of system capability.

B.01.05 Release Note

What's New In Agilent Headspace Drivers Version B.01.05

Compatibility with Agilent Data System

Supported Agilent Headspace Hardware

Compatibility with Operating System

What's' New in Agilent Headspace Drivers Version B.01.05

Resolved issues reported by customers. Upgrades will be covered via this release note if customers have experienced these issues.

Issues Resolved in B.01.05:

| KPR | Title |
|-----|--|
| 123 | G1888 Method Assist 'Use existing method timing' does not retain the GC Cycle Time. |
| 124 | Headspace does not go to Sleep as scheduled. |
| 125 | Error 1330 occurs during Headspace Driver installation. |
| 129 | 7697 Detailed Status does not number the extraction/injections when run in Concentrated Extraction Mode. |
| 130 | 7697 Detailed Status does not number the injections when run with MHE mode. |

Compatibility with Agilent Data Systems

Headspace software, revision B.01.05 was tested on /supported with,

| Software | Revisions |
|-------------------------|-------------------|
| GC ChemStation | B.04.03 SP1 |
| OpenLAB CDS ChemStation | C.01.01-C.01.05 |
| OpenLAB CDS EZChrom | A.04.02 – A.04.05 |
| GC MassHunter | B.07.00 |
| GCMS ChemStation | E.02.02 SP2 |

Supported Agilent Headspace Hardware

Headspace software, revision B.01.05 was tested on /supported with,

7697A Headspace Hardware

| Model | Description | FW Revision | ICF Ready Package | Communication |
|--------|--------------------------------------|-------------|-------------------|---------------|
| G4556A | 7697A Headspace Sampler, 12 vial | A.01.05.1 | A.02.02 | TCP/IP |
| G4557A | 7697A Headspace Sampler, 111 vial | A.01.05.1 | A.02.02 | TCP/IP |

G1888A Headspace Hardware

| Model | Description | FW Revision | ICF Ready Package | Communication |
|--------|-------------|-------------|-------------------|---------------|
| G1888A | Headspace | A.01.10 | A.02.02 | TCP/IP |

Compatibility with Operating Systems

| Components List | System Type | Revision | Language |
|-------------------------|-------------|----------|---|
| Windows 8 Enterprise | 64-bit | | English US, Japanese, Chinese, Brazilian Portuguese |
| Windows 8 Enterprise | 32-bit | | English US, Japanese, Chinese, Brazilian Portuguese |
| Windows Server 2008 | 64-bit | R2 | English US, Japanese, Chinese, Brazilian Portuguese |
| Windows Server 2003 | 32-bit | SP3 | English US, Japanese, Chinese, Brazilian Portuguese |
| Windows 7 Enterprise | 32 Bit | SP1 | English US, Japanese, Chinese, Brazilian Portuguese |
| Windows 7 Enterprise | 64 Bit | SP1 | English US, Japanese, Chinese, Brazilian Portuguese |
| Windows XP Professional | 32 Bit | SP3 | English US, Japanese, Chinese, Brazilian Portuguese |