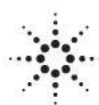




Agilent OpenLAB Data Store

Version A.02.02

**Release Notes and
History**



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2014

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

M8620-90114

Edition

October 2014

Agilent Technologies, Inc.

Software Revision

This guide is valid for Agilent OpenLAB Data Store version A.02.02.

Microsoft ® is a U.S. registered trademark of Microsoft Corporation.

This product may be used as a component of an in vitro diagnostic system if the system is registered with the appropriate authorities and complies with the relevant regulations. Otherwise, it is intended only for general laboratory use.

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

If software is for use in the performance of a U.S. Government prime contract or subcontract, Software is delivered and licensed as "Commercial computer software" as defined in DFAR 252.227-7014 (June 1995), or as a "commercial item" as defined in FAR 2.101(a) or as "Restricted computer software" as defined in FAR 52.227-19 (June 1987) or any equivalent agency regulation or contract clause. Use, duplication or disclosure of Software is subject to Agilent Technologies' standard commercial license terms, and non-DOD Departments and Agencies of the U.S. Government will receive no greater than Restricted Rights as defined in FAR 52.227-19(c)(1-2) (June 1987). U.S. Government users will receive no greater than Limited Rights as defined in FAR 52.227-14 (June 1987) or DFAR 252.227-7015 (b)(2) (November 1995), as applicable in any technical data.

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.

Introduction

This document describes the functions, features and changes introduced with this and previous releases of the Agilent OpenLAB Data Store product.

References to product documentation regarding known issues and workarounds are also provided.

For our regulated customers:

When any change is made to Agilent software, the validation status of the software needs to be re-established by the user/customer. Whenever software is changed, a validation analysis should be conducted not just for the validation of an individual change, but also to determine the extent and impact of that change on the entire software system. Agilent provides a document that can help with the impact analysis.

Version A.02.02

New Operating System Support

- Microsoft Windows Server 2012 R2

New Databases Support

- PostgreSQL 9.2
- Microsoft SQL Server 2012 Service Pack 2

New Features

- FTP Support
Generic FTP client tools like FileZilla can be used to upload files to the Data Store repository or to download files. The user's permissions and privileges configured on the system are also applicable for the FTP activities.
The FTP protocol is enabled by default, the procedure to enable or disable FTP is documented in the OpenLAB Data Store Administrator Guide.
- Support of HTTPS and FTPS

Version A.02.01

The OpenLab Data Store A.02.01 server can be installed independent of other OpenLAB software through a dedicated installer program.

The software can be installed on a PostgreSQL database engine rev. 9.0 in addition to Microsoft SQL Server 2008 R2 SP2.

The server is automatically activated upon installation with internal authentication. A default user called 'admin' with the password 'openlab' is created.

The software supports up to 30 instruments of mixed OpenLAB CDS WorkStations (ChemStation or EZChrom) and ICP-MS MassHunter WorkStations on the same server.

Lab Applications introduced with OpenLAB Data Store A.01.02 is not supported with this release.

Version A.01.03

OpenLAB Data Store A.01.03 “*OpenLAB Data Store for MS*” ONLY supports the Agilent ICP-MS MassHunter software G7201B rev. B.01.03.

This release introduces the support of the PostgreSQL database engine rev. 9.0 in addition to Microsoft SQL Server 2008 R2 SP2.

Lab Applications introduced with OpenLAB Data Store A.01.02 are not supported for OpenLAB Data Store for MS.

Version A.01.02

With OpenLab Data Store A.01.02, the LAB Applications have been introduced to accommodate common laboratory workflows and include two applications:

Sample Submission

- Request an analysis - Prepare a chromatography sequence and assign it to the analyst through Data Store.
- Accept and complete the analysis - The analyst receives the sequence, runs it on an instrument using an Agilent CDS system, and stores the results into Data Store.
- Review results - Import the results, review, approve or reject them in Data Store.

Lab Journal

- Manage lab assets, such as instruments and columns.
- Capture important lab events such as instrument service, replacements, calibrations and etc.
- Review the operations that occurred on a specific instrument or column.
- Display simple usage statistics for instruments or columns.

In addition, the Data Store user interface has been redesigned to support the new applications.

Version A.01.01

OpenLAB Data Store A.01.01 Supported with CDS

OpenLAB Data Store is a central repository where you can safely archive, store, and share electronic files.

OpenLAB CDS ChemStation Edition and Shared Services support the use of OpenLAB Data Store as the storage location for your scientific data. Use of Data Store storage is supported in networked and distributed configurations of OpenLAB CDS and provides several benefits for your system:

- Central Storage with security for small to medium sized labs
- Handles localized Chinese/Japanese content
- Web based data access for search, share, and review
- 21 CFR 11 Compliant e-signatures and audit trail

Known Issues

For known issues and workarounds in the OpenLAB Data Store software at the time of release see the Software Status Bulletin file named *M86xxAA-SSB.html* located on Disk 1 of the OpenLAB Data Store Installation media.

For known issues that have been resolved and/or closed with a workaround, see the Software Release Bulletin file named *M86xxAA-SRB.html* located on Disk 1 of the OpenLAB Data Store Installation media.

You may also visit www.Agilent.com for up to date issue information for all Agilent software products:

<http://www.chem.agilent.com/en-US/Support/Downloads/Patches/Pages/default.aspx>