



Agilent Dissolution Workflow Manager (DWM)
for OpenLab

System Requirements

Notices

Manual Part Number

Document No: D0125049 Rev. A.00
Edition 02/2025

Copyright

© Agilent Technologies, Inc. 2025

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.

Agilent Technologies, Inc.
5301 Stevens Creek Blvd.
Santa Clara, CA 95051
USA

Software Revision

This guide is valid for version 1.0 of OpenLab DWM.

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.

Contents

1	Introduction	4
	Terms and Abbreviations	5
	Version History	5
2	OpenLab DWM Topologies	6
	OpenLab DWM Workstation Plus	6
	OpenLab DWM Client/Server	7
3	PC Requirements	9
4	Configuration Capacity	10
5	Instrument Requirements	11
	Supported Instruments	11
6	Licensing	12
7	Known Issues and Limitations	13
8	Troubleshooting	14
	Initial Troubleshooting Steps	14
	Using Online Help	14
	Support	15

Introduction

This document details the additional minimum hardware and software requirements that need to be met to run DWM for OpenLab. It is valid for Workstation Plus, client, or analytical Instrument Controller (AIC) components. It also provides information on supported Agilent Instrumentation.

A full installation of OpenLab CDS v2.8 update 5 is required prior to installation of Dissolution Workflow Manager (DWM) for OpenLab, see [documents for Installation and Configuration of OpenLab CDS](#) for details.

Terms and Abbreviations

Table 1 Terms and Abbreviations

Term	Description
OpenLab Control Panel	Control panel for Agilent OpenLab Software
Microsoft Control Panel	Part of the Microsoft Windows operating system
CDS	Chromatography Data System
AIC	Agilent Instrument Controller – part of a distributed Client/Server system
DWM	Dissolution Workflow Manager
FW	Firmware

Version History

Table 2 Software Version History

Software Version release date	Version revision number	Description
February 2025	1.0	Initial Release

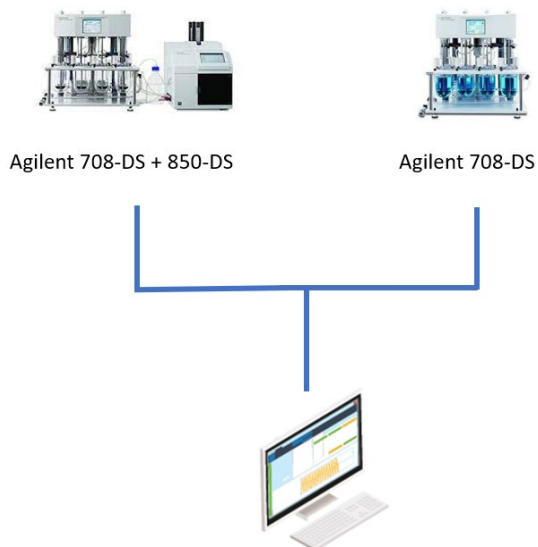
2

Dissolution Workflow Manager (DWM) for OpenLab Topologies

Dissolution Workflow Manager (DWM) for OpenLab is available in different topologies. They differ in the location and availability of OpenLab software components (shared services, secure storage, clients). For details on the topologies refer to the respective installation guide or check with your Agilent representative.

Dissolution Workflow Manager (DWM) for OpenLab Workstation Plus

Dissolution Workflow Manager (DWM) for OpenLab Workstation Plus installations include all required components on the same PC. Workstation Plus installations include content management (provided by the local Secure Storage component).



NOTE

Standard OpenLab CDS Workstation installation are not compatible with OpenLab CDS.

Dissolution Workflow Manager (DWM) for OpenLabClient/Server

Dissolution Workflow Manager (DWM) for OpenLab installs components onto the client PCs, Analytical Instrument Controllers (AICs) and Servers.

Consult with your Agilent support representative to decide which topology and requirements are appropriate for your environment.



3 PC Requirements

Dissolution Workflow Manager (DWM) for OpenLab requires one RS-232 port per instrument module to be connected to the Workstation Plus PC or Analytical Instrument Controller (AIC). Consult with your agilent representative to discuss USB to RS-232 module options available should there be insufficient RS-232 ports available. For all other PC requirements, consults the CDS v2.8 System requirements guide.

4 Configuration Capacity

Instrument: An analytical system, made up of one or more modules that are configured together to achieve the desired analytical functionality. Also known as a system, i.e. an LC/MSD system.

Points: The relative level of load (load value) that an instrument or module puts on a physical or virtual system, aggregating CPU and RAM (memory) consumption. Different types of detectors require more processing power and memory than others, so they have been assigned higher “point” values.

Concurrent sessions: The number of simultaneous connections to the system. This includes user connections and instrument connections. For OpenLab CDS, use points as the measure of instrumentation as this reflects both number of instruments and the amount of data generated by each instrument system. See below for examples of number of points per instrument type.

Instrument licenses: Connection Licenses refer to the licenses required in the software in order to run a specific instrument configuration. Note: Points do not equal instrument connection licenses.

Dissolution Workflow Manager (DWM) for OpenLab instrument control is operated via Dissolution Workflow Manager (DWM) for OpenLab counted licenses. Therefore, there is no limit to the number of instruments configured on Dissolution Workflow Manager (DWM) for OpenLab. Only the number of instruments connected at a given time. A license is returned to the License Pool following instrument connection closure.

An individual Workstation Plus or AIC is capable of processing 4 connection points or 2 instrument connections concurrently.

Instrument Configuration	Number of Licenses Consumed	Instrument Points Consumed
708-DS Standalone	1	1
708-DS / 850-DS System	1	2

5 Instrument Requirements

Supported Instruments

Dissolution Workflow Manager (DWM) for OpenLab supports the following dissolution instruments:

Instrument Module	Module Type	System FW
708-DS	Dissolution Bath	3.01 or greater
850-DS	Autosampler	4.01 or greater
850-DS w/ filter changer	Autosampler	4.01 or greater

To obtain firmware updates for the relevant instruments, contact your Agilent support representative.

6

Licensing

Dissolution Workflow Manager (DWM) for OpenLab operates using dedicated Dissolution Workflow Manager License, this can be generated via SubscribeNet.

Dissolution Workflow Manager (DWM) for OpenLabinstrument control is operated via Dissolution Workflow Manager (DWM) for OpenLabcounted licenses. Therefore, there is no limit to the number of instruments configured on Dissolution Workflow Manager (DWM) for OpenLab. Only the number of instruments that can be connected at a given time. A license is returned to the License Pool following instrument connection closure.



7

Known Issues and Limitations

Consult the Dissolution Workflow Manager (DWM) for OpenLab SSB for a list of all current known issues and limitations.

This can be found on the installation USB and via the Agilent Dissolution Community page.

Initial Troubleshooting Steps

- ✓ Double-check that all instrument components are connected and turned on.
- ✓ Verify that the instruments and modules you have connected are supported (see "Supported Instruments and Modules" on page 11).
- ✓ Verify that the firmware revision meets the minimum requirements for the specified instrument.
- ✓ Double-check the LAN/IP address settings as documented in this guide.
- ✓ Check that the driver is correctly installed and reports the correct version by checking what is listed in Windows **Programs and Features** (or equivalent) or running the Software Verification tool.
- ✓ Check that all relevant updates and hotfixes for the driver have been applied.
- ✓ If you installed the software on an OpenLab CDS Distributed System You have properly set up all additional DCOM settings and Allow permissions on the AIC. Refer to Table 8 on page Fehler! Textmarke nicht definiert. and Table 9 on page Fehler! Textmarke nicht definiert..

Using Online Help

Online help can be accessed using the ? icon while the user interface is open within the Agilent OpenLab Software.

Support

For support with this software, contact via submitting a ticket through the support portal:

<https://www.agilent-labinformatics.com/>

In This Book

This installation guide is designed to help users install and configure instruments for use with Agilent Dissolution Workflow Manager (DWM) for OpenLab software.

The manual describes the following:

- Introduction
- Requirements
- Installation
- Configuration
- Uninstall the Drivers
- Troubleshooting

www.agilent.com

© Agilent Technologies, Inc. 2025

Edition 02/2025

Document No : D0125049 Rev :A.00

