



ADFExport for OpenLab



Notices

Document Information

Document No: D0006043
Edition 09/2021

Copyright

© Agilent Technologies, Inc. 2020-2021

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 1.3 of ADFExport for OpenLab.

Product No: M8480AA

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.

In this Guide ...

This guide describes the requirements and installation of ADFExport for OpenLab rev. 1.3.

ADFExport for OpenLab was designed to work with two Agilent Chromatography Data Systems: OpenLab CDS and OpenLab CDS ChemStation Edition. Some features or behaviors of the ADFExport plugin differ depending on your system. This document includes details on these differences.

Table 1 Terms and abbreviations

Term	Description
CDS	Chromatography Data System
ChemStation	OpenLab CDS ChemStation Edition
Content Management	Data storage component provided as part of the OpenLab system
OpenLab CDS	Version 2.x of OpenLab CDS
Control Panel	Control Panel for Agilent OpenLab software
Microsoft Control Panel	Part of the Microsoft Windows operating system
AIC	Agilent's Analytical Instrument Controller

Contents

1	Introduction	6
	ADF File Content	6
2	Licensing	7
	Available Products	7
	Get a License	7
	Install Your License	8
3	ChemStation	9
	Requirements	9
	Installation	10
	Get Installation Package	10
	Prepare Network Drives	10
	Install ADFExport for OpenLab	10
	Software Verification Tool	11
	Verify Network Access Privileges	11
	Upgrade ADFExport for OpenLab	12
	Backup customized Macros	12
	Upgrade Instructions	12
	Software Migration	14
	Re-register Macro in the Run Time Checklist	14
	Move Data to the New System	14
	Data Migration	15
	Uninstallation	16
4	OpenLab CDS	17
	Requirements	17
	Installation	18
	Get Installation Package	18
	Prepare Network Drives	18
	Install ADFExport for OpenLab	18

Contents

Software Verification Tool	19
Verify Network Access Privileges	19
Upgrade ADFExport for OpenLab	20
Uninstallation	20
Appendix: License Terms	21

Introduction

ADFExport for OpenLab is an Add-on that offers the possibility to export results to the Allotrope Data Format (ADF). A detailed description of the Allotrope Data Format and its structure is available on the Allotrope Foundation web site (<https://www.allotrope.org>).

ADF File Content

In the current revision of the ADFExport Add-on, the following content is written into the ADF file (for detailed information see **ADFExport_ADFContentDescription.pdf** available in the download package in SubscribeNet).

- ADF data description

The data description contains an initial set of metadata about the single run or the entire sequence. The metadata includes information about the run, such as the system (instrument, including module and column information), the submitter, the sample, the injection, and the sequence. Furthermore, information on chromatograms and UV spectra as well as acquisition and processing method names and peak results are stored.

- ADF data cube

The data cube contains exported raw data (chromatograms and spectra).

- ADF data package

The data package contains a subfolder with the original OpenLab CDS or ChemStation files.

For more information about the ADF file content, see the ADFExport for OpenLab online help.

2

Licensing

Every chromatography data system that runs with an ADFExport Add-on requires a license. Each time the ADFExport Add-on is used, it checks the licensing server for an available license and uses it during its operation. If the ADFExport Add-on is stopped or fails, the license is released again.

The number of licenses depends on the number of ADFExport instances you want to use at the same time.

For information on the License terms, see “[Appendix: License Terms](#)” on page 21.

Available Products

Table 2 Available Products for ADFExport for OpenLab

Product number	Product name	Description
M8480AA	ADFExport for OpenLab rev. 1.3	Supports one software instance (1 instrument)

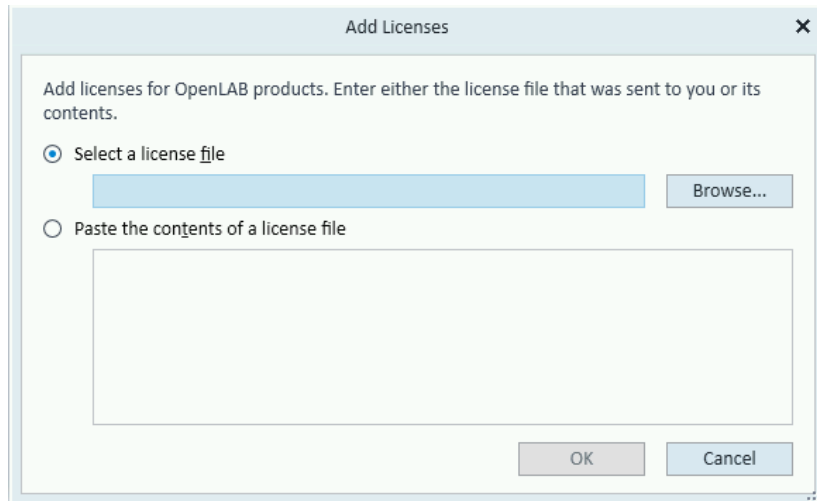
Get a License

Log in to SubscribeNet to generate a license. During this process you will have to enter the MAC address of your license server. For client/server systems, this is the server. Download your license file.

For more details, please refer to the *OpenLab CDS Clients and Instrument Controllers* guide (CDS_ClientAIC.pdf) or *OpenLab CDS ChemStation Edition Networked and Distributed System Installation* guide (CDS_CS_Install_NwWS-DS.pdf).

Install Your License

- 1 In the OpenLab Control Panel under **Administration**, select **Licenses** and click **Add License**.



The screenshot shows a dialog box titled "Add Licenses" with a close button (X) in the top right corner. The main text reads: "Add licenses for OpenLAB products. Enter either the license file that was sent to you or its contents." There are two radio button options. The first option, "Select a license file", is selected and has a text input field followed by a "Browse..." button. The second option, "Paste the contents of a license file", is unselected and has a large text area below it. At the bottom right of the dialog are "OK" and "Cancel" buttons.

- 2 Choose to install the license by:
 - Using the license file option to browse to and open the license file (.lic) saved from the license generation process in SubscribeNet.
 - Selecting the License Text option and copying the license text from a text file received into the provided field.
- 3 Click **OK**.

The Administration interface in the Control Panel will now display the status of installed licenses.



3

ChemStation

Requirements

ADExport for OpenLab, rev. 1.3, is compatible with ChemStation C.01.10 and C.01.09 Update 2.

Signals from Agilent ChemStation LC (SFC or HPLC) instrument runs acquired with DAD, VWD or MWD can be exported.

ADExport does not support to export ChemStation data acquired from hybrid SFC/HPLC systems nor from third-party vendor LC, LC/MS, GC/MS, nor GC signals.

The product is not localized and can only be used on English versions of ChemStation.

Installation

Get Installation Package

Download the **ADFExport for OpenLab** installation package from SubscribeNet, and extract the files to a local folder or to a network drive.

Prepare Network Drives

If you will use the ADFExport installation package from a network drive, you must prepare the network drive to make it accessible by the wizard. Without this preparation, Windows security prevents the wizard from accessing those drives.

- 1 Map the drive to a letter.

For example, map the drive as **Z:** using the shared path "\\<machine-name>\OpenLabCDS".

This maps the drive for the logged-in user.

- 2 Open the command prompt in elevated mode (run as administrator), and map the drive using the net use command.

For example,

```
net use Z: "\\<machine-name>\OpenLabCDS"
```

This maps the drive for the local administrator account. The mapped drive is now visible to both logged-in user and administrator, and can be selected in the wizard.

Install ADFExport for OpenLab

Install the software on the PC on which your instruments are configured, that is, the Workstation or Networked Workstation. In Distributed Systems, install it on the Analytical Instrument Controller (AIC).

- 1 Go to **Start > All programs > Agilent Technologies > OpenLab Additional Software and Drivers**.
- 2 On the **Planning** page, select the **Install** option.
- 3 On the **Add-on Selection** page, select the **Agilent.OpenLab.ChemStation.ADFExport.xml** file of your installation package.
- 4 Follow the wizard to install the required software.
- 5 Shut down and restart your ChemStation instances to make the add-on available.

Software Verification Tool

The Software Verification Tool (SVT) runs automatically after installing ADFExport for OpenLab via **OpenLab Additional Software and Drivers**.

The SVT report informs you about successful integration into the existing ChemStation system. The single files installed by the ADFExport Add-on are not listed in the SVT report.

Verify Network Access Privileges

If ADF files shall be exported to a network share, users need sufficient privileges to access the share. In a Distributed System, these privileges must be granted to the AIC Instrument user.

Upgrade ADFExport for OpenLab

Backup customized Macros

If you made any changes to the original **CSADFExportSingleSampleRun.mac** file provided with ADFExport for OpenLab rev. 1.0, or used this file for any other purposes than the default export: Please make sure to make a safety copy of the macro files. The macro file **CSADFExportSingleSampleRun.mac** will be removed during the upgrade.

Upgrade Instructions

- 1 If you run ChemStation C.01.09 with ADFExport for OpenLab rev. 1.0; Upgrade ChemStation to version C.01.09 Update 2 or higher as described in the ChemStation manuals (upgrading ADFExport rev. 1.0 to 1.3 is not supported on lower ChemStation revisions).
- 2 Go to **Start > All programs > Agilent Technologies> OpenLab Additional Software and Drivers**.
- 3 On the **Add-on Selection** page, select the **Agilent.OpenLab.ChemStation.ADFExport.xml** file of your installation package.
- 4 A confirmation dialog opens. Confirm that you want to upgrade the older version of the ADFExport Add-on.
- 5 Follow the wizard.

Default Output Path

If you upgrade ADFExport for OpenLab from rev. 1.0, the default output path for automated ADF exports changes.

In rev. 1.0, it was **C:\Users\Public\Documents\ChemStation\<instrument>\data**.

After successful upgrade, it is set to **C:\Users\Public\Documents\ChemStation\<instrument>\CSADFExport**.

If you upgrade from rev. 1.2, the default output path does not change.

Default Log File Location

If you upgrade ADFExport for OpenLab from rev. 1.0, the default path for log files created by exports from the Windows Command Prompt changes.

In rev. 1.0, it was

\Users\<user name>\Documents\CSADFExportLogs.

After successful upgrade, it is set to

%ProgramData%\Agilent\LogFiles\CSADFExportLogs.

If you upgrade from rev. 1.2, the log file location does not change.

To Adjust the Output Path per Instrument (ChemStation C.01.10)

- 1 Open the Agilent OpenLab Control Panel.
- 2 Select the **Instruments** tab.
- 3 Navigate to the relevant instrument, and click the **Configure Instrument** icon or right click the instrument name and select **Configure Instrument**.

The **Configure Instrument** dialog box is displayed.

- 4 In the bottom-right corner, select **Additional configuration > Paths...**

The **Configure Path Variables** dialog box is shown.

- 5 For the **ADF Output Path**, enter the required path.
- 6 Confirm your changes.

The changed output path will automatically be used next time when you run the **Export to ADF** Macro/Command.

To Adjust the Output Path per Instrument (ChemStation C.01.09 Update 2 or higher)

- 1 Open a ChemStation instance for the relevant instrument.
- 2 If the command line is not already visible, select **View > command line** in the **Method and Run Control** view to show it.
- 3 Enter the command `SetADFOutputPath "<newPath>"`.
- 4 For example: `SetADFOutputPath "C:\Users\Public\Documents\ChemStation\1\my_CSADFExport\"`
- 5 The default output path is changed in the ChemStation.ini file for the current instrument

Software Migration

Re-register Macro in the Run Time Checklist

This procedure is required if the following applies:

- You upgrade ADFExport for OpenLab from rev. 1.0.
- You used the post-run command/macro *as provided by ADFExport*, and you continue using the relevant ChemStation methods.

Update the ChemStation methods as follows:

- 1 In ChemStation, load the relevant method.
- 2 Open the Run Time Checklist.
- 3 Under **Post-Run Command / Macro**, delete the command name field.
- 4 Clear the **Post-Run Command / Macro** check box.
- 5 To close the dialog, click **OK**.
- 6 Open the Run Time Checklist again.
- 7 Select the **Post-Run Command / Macro** check box.
- 8 From the drop-down list for post-run commands, select **Export to ADF**.
- 9 Confirm your changes.
- 10 Save the updated method.

Move Data to the New System

If you upgrade ADFExport for OpenLab from rev. 1.0: To move your data from the old to the new system, transfer the data to the output path for automated ADF exports. For location details, see:

- ["Default Output Path"](#) on page 12
- ["To Adjust the Output Path per Instrument \(ChemStation C.01.10\)"](#)
- ["](#) on page 13

Data Migration

If you have generated an ADF file with rev 1.2 or lower of ADFExport forOpenLab, you can update its content to the current format. To update an ADF file, use the Windows command prompt.

Prerequisites: Ensure that the ADF file name does not include any of the following characters - they are not supported anymore: ` ^ % # °`

- 1 From the Windows Start menu, select Windows System > Command Prompt.

The Windows Command Prompt window opens.

- 2 Use the *cd* (change directory) command to change to the **CSADFExport** subfolder in the ChemStation installation folder.

By default, the path is **C:\Program Files (x86)\Agilent Technologies\CSADFExport**. To go there, type *cd C:\Program Files (x86)\Agilent Technologies\CSADFExport* and press **[Return]**. If your ChemStation installation folder is different from the default, replace the **\Program Files (x86)\Agilent Technologies** with the appropriate path definition.

The **CSADFExport** folder contains the executable file, **Agilent.OpenLab.CSADFExport.exe**.

- 3 To migrate the ADF file, type *Agilent.OpenLab.CSADFExport.exe -m basic -i "<input path>" -o "output path"* at the command prompt.

-i specifies the input path. It is the full path to the existing ADF file, including the file name itself.

-m specifies the export mode. For the current release, the only valid argument is basic.

-o specifies the output path. If the specified path does not exist or **-o** is not specified, the ADF file is stored in **C:\Users\<user name>\Documents\CSADFExport**. To export files to a network share, enter the output path using UNC syntax (**\\hostname\share\path**).

Uninstallation

To uninstall ADFExport for OpenLab:

- 1 Go to **Start > All programs > Agilent Technologies> OpenLab Additional Software and Drivers**.
- 2 On the **Installed Features** page, select **Agilent ADFExport for OpenLab**.
The uninstallation wizard opens.
- 3 On the **Planning** page, select the **Uninstall** option.
- 4 Follow the wizard to uninstall the plug-in.

4

OpenLab CDS

Requirements

ADFExport for OpenLab, rev. 1.3, is compatible with OpenLab CDS 2.4, 2.5, and 2.6.

Signals from Agilent OpenLab CDS LC (SFC or HPLC) instrument runs acquired with DAD, VWD or MWD can be exported.

ADFExport does not support to export CDS data acquired from hybrid SFC/HPLC systems nor from third-party vendor LC, LC/MS, GC/MS, nor GC signals.

The product can be used on Chinese, Japanese, Portuguese, or Russian versions of OpenLab CDS.

Installation

Get Installation Package

Download the **ADExport for OpenLab** installation package from SubscribeNet and extract the files to a local folder or to a network drive.

Prepare Network Drives

If you will use the ADExport installation package from a network drive, you must prepare the network drive to make it accessible by the wizard. Without this preparation, Windows security prevents the wizard from accessing those drives.

- 1 Map the drive to a letter.

For example, map the drive as **Z:** using the shared path "\\<machine-name>\OpenLabCDS".

This maps the drive for the logged-in user.

- 2 Open the command prompt in elevated mode (run as administrator), and map the drive using the net use command.

For example,

```
net use Z: "\\<machine-name>\OpenLabCDS"
```

This maps the drive for the local administrator account. The mapped drive is now visible to both logged-in user and administrator, and can be selected in the wizard.

Install ADExport for OpenLab

Install the software on the PC on which your instruments are configured. In Client/Server systems, install it on the Analytical Instrument Controller (AIC) and on all relevant clients.

- 1 Run **Agilent.OpenLab.CDS.ADExport.msi** as an administrator.
- 2 Follow the wizard.
- 3 Install the required license (see "Install Your License" on page 8).

Software Verification Tool

To qualify the ADFExport for OpenLab installation, launch the Agilent Software Verification Tool (from Agilent Technologies/Software Verification Tool).

The SVT report informs you about successful integration into the existing OpenLab CDS system. The SVT report includes the files installed by the ADFExport Add-on.

Verify Network Access Privileges

If ADF files shall be exported to a network share, users need sufficient privileges to access the share. In a Distributed System, these privileges must be granted to the AIC Instrument user.

Upgrade ADExport for OpenLab

Simply run the installation wizard as described under “Install ADExport for OpenLab” on page 18. You do not need to uninstall ADExport for OpenLab.

If you are using ADExport for OpenLab and plan to upgrade OpenLab CDS:

- 1 Uninstall the ADExport Add-on.
- 2 Upgrade OpenLab CDS.
- 3 Install the current revision of ADExport for OpenLab.

NOTE

ADExport for OpenLab from rev. 1.1 is not supported anymore with OpenLab CDS revisions higher than 2.4, and rev.1.2 is not supported with OpenLab CDS revisions higher than 2.5. If you upgrade OpenLab CDS, you must also upgrade ADExport for OpenLab from rev. 1.1 or 1.2 to rev. 1.3. To upgrade please follow step 1 to 3.

Uninstallation

- 1 Log in as an administrator.
- 2 In the Windows settings, open **Apps > Apps & features**.
- 3 Select **ADExport for OpenLab**, and confirm to uninstall it.
- 4 Reboot.



Appendix: License Terms

AGILENT TECHNOLOGIES, INC. SOFTWARE LICENSE TERMS FOR Agilent OpenLAB Software

For AGILENT TECHNOLOGIES, INC. SOFTWARE LICENSE TERMS FOR Agilent OpenLab Software read OpenLab_LicenseTerms.pdf provided with installation.

Allotrope Foundation ADF license information

ADFExport software makes use of and incorporates Allotrope Foundation ADF framework. Please read Allotrope_ADFLicenseInformation.pdf provided with installation.

commandlineparser/commandline license information

ADFExport software makes use of and incorporates commandlineparser/commandline. Please read below License and copyright notice.

The MIT License (MIT)

Copyright (c) 2005 - 2015 Giacomo Stelluti Scala & Contributors

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

In This Book

This manual describes the ADFExport for OpenLab CDS rev. 1.3:

- OpenLab CDS ChemStation Edition and OpenLab CDS
- Licensing
- Requirements
- Installation
- Upgrade
- Migration
- Uninstallation

www.agilent.com

© Agilent Technologies, Inc. 2021

Edition 09/2021
Document No: D0006043

