

Migration From Agilent Alliance Driver To Driver Central Alliance Driver

Purpose: This document is for customers who are currently using the Agilent Alliance driver with OpenLAB CDS EZChrom Edition and will need to move to the Driver Central Alliance driver. It describes the procedure for migrating from the Agilent Alliance driver to the Driver Central Alliance driver.

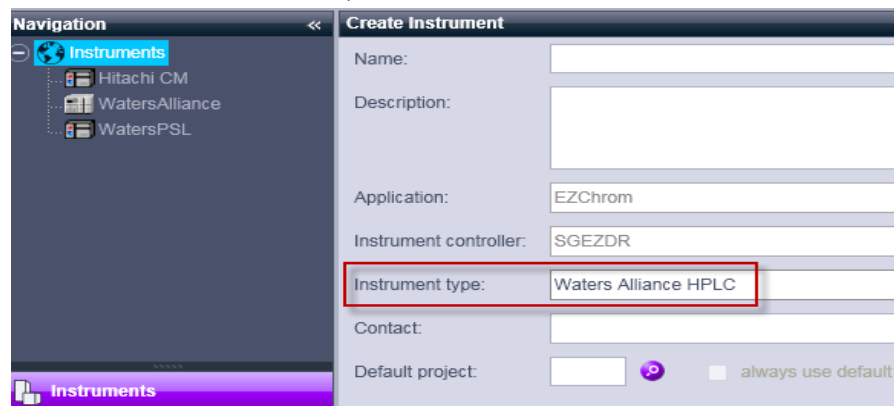
Use of the Agilent Alliance Driver: The Agilent Alliance driver should be used if you want to reprocess data acquired with the Agilent Alliance driver. Data acquired with the Agilent Alliance driver can be viewed in the Driver Central Alliance driver as long as there is no re-processing and there is no need to view your method. When you open the data a message will be displayed indicating that you are loading a method that does not match the instrument configuration. Select “No” when prompted to save the Method. All new Acquisitions need to be performed with the Driver Central Alliance driver. Data acquired with the Driver Central Alliance driver cannot be reprocessed or viewed using the Agilent Alliance driver.

Having both the Agilent Alliance Driver and Driver Central Alliance Driver on the same machine:

Both the Agilent Alliance driver and the Driver Central Alliance driver can co-exist on the same machine. Only the Driver Central Alliance driver will be used with online instruments for data acquisition.

Migration From Agilent Alliance Driver to Driver Central Alliance Driver:

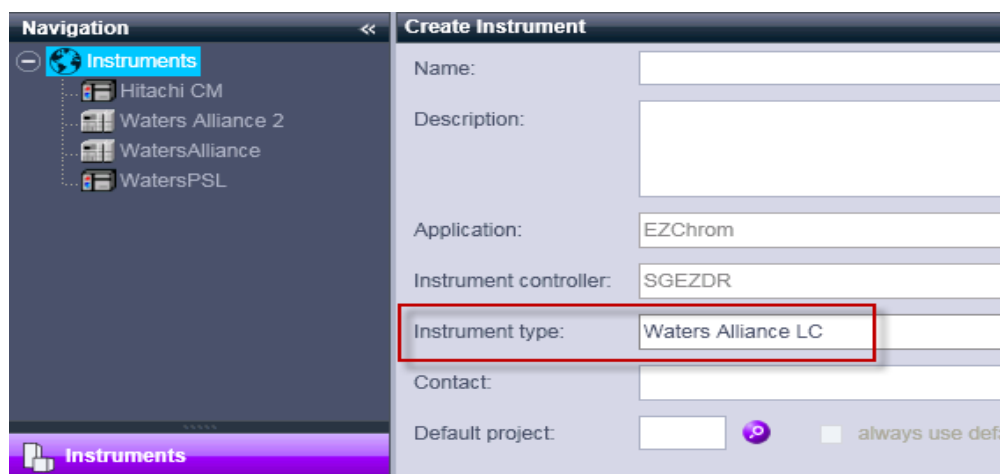
1. Install the Driver Central Alliance driver using the provided instructions.
2. Create and configure a new instrument with the Driver Central Alliance driver. Select “Waters Alliance HPLC” to create an instrument with the Driver Central Alliance driver. This will be the instrument used for data acquisition.



3. Create and configure a new instrument with the Agilent Alliance driver. Select “Waters Alliance LC” to create an instrument with the Agilent Alliance Driver. When you configure the instrument you do not need to make an instrument connection since all work with this driver will be performed using an offline session. To configure the Agilent Alliance driver do the following:
 - a. Click “Configure Instrument”.
 - b. Move the desired modules from “available modules” to “configured modules”.

- c. Open the modules.
- d. Click OK without editing or making any changes.

Now you will have instrument that will be used only for offline sessions to perform data analysis.



4. New Methods will need to be created with the Driver Central Alliance driver. All Agilent Alliance driver methods will need to be recreated in the Driver Central Alliance driver. Use the procedure below to migrate your Agilent Alliance Methods to the Driver Central Alliance Methods.
 - a. Open an Agilent Alliance instrument and a Driver Central Alliance instrument offline.
 - b. Open your Agilent Alliance Method.
 - c. Create a new Driver Central Alliance method using the Agilent Alliance method opened in step b. New Method creation is a manual process.

Differences Between the Agilent Alliance Driver and Driver Central Alliance Driver:

The table below describes the differences between the Agilent Alliance driver and the Driver Central Alliance driver.

	Agilent Alliance Driver	Driver Central Alliance Driver
Data Collection Mode for 2996 DAD Detector	3D Only	2D and 3D
Spectra Resolution for PDA	2nm, 3nm, 4nm...	1.2nm, 2.4nm, 3.6nm
Signal Unit	AU	mAU

Temperature Control	Temperature range (for readiness) is hardcoded to 5°C.	Temperature range (for readiness) is defined in Method.
Threshold program on 2487 UV Detector	Not Supported	Supported
Intelligent Reporting	Not Supported	Supported