

Agilent PC Bundle for Chemstation Edition

PC Bundle Set-up with Microsoft Windows 10

First time boot-up of the Agilent bundle PC

- 1 When you start this PC the first time, an internal script is activated. The user is prompted to select the country/region, preferred language, keyboard layout and the time zone, user name and password.

NOTE

The language must be set to be one of the Agilent supported Operating System languages English (United States), Chinese or Japanese

- 2 As soon as the Microsoft Windows 10 related configuration is done, and the desktop is visible **restart the Bundle PC to complete the Windows setup before doing the next steps!** (This may take several minutes)
- 3 The **PC name** will now be DESKTOP-XXXXXXX (where "XXXXXXX" is a randomly generated sequence of letters and numbers). If a different PC name is required change it now. If a PC name change is required after ChemStation got installed, please check the installation manual for further instructions.
- 4 Connect the PC to a given domain if desired.
- 5 **Ensure all Windows Updates have completed before installing ChemStation.**

- 6 Click the **OpenLabCDSCS** icon  on the desktop to run the install script.

Don't close the black CMD window during the installation process.

After the installation the Software Verification Report will show up. This indicates that the installation is done.

NOTE

In case the CMD window is still open it can be closed now.

- 7 Open the OpenLab Control Panel  and install your license. The OpenLab CDS ChemStation Edition is now ready to use.

Help and Support

Using Online Help

Online Help can be accessed using via **F1** within any open user interface interacting with Agilent OpenLab Software.

Getting Support

Please contact your local Agilent Technologies Customer Care Center to obtain support. For details please visit the following web page: <http://www.agilent.com/chem/contactus>.

Agilent Technologies also provides electronic support through the World Wide Web via <http://www.agilent.com/chem>.

In any communication with the Software support team regarding a problem with this OpenLab CDS bundle, please clearly state the following:

- Your name, address, e-mail address and telephone number.
- Your OpenLab CDS product name and version number. Both are listed in Help and About.
- A description of the problem, and what you were trying to do when the problem occurred.

Re-Image and Installation Preparation Procedure

This section refers to the Emergency Recovery Media provided with the **HP Z2-G4 Workstation Bundle PC for OpenLab CDS ChemStation** from Agilent.

The system will be set to the following configuration:

Edition: Windows 10 Pro
Version: 1803
OS Build: 17134.1
System type: 64-bit operating system

CAUTION Irreversible loss of data will occur when re-imaging the computer!

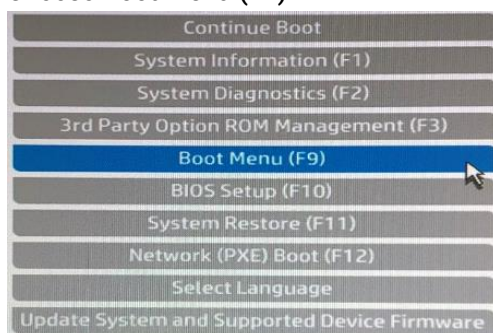
STEPS FOR RE-IMAGING THE PC

NOTE

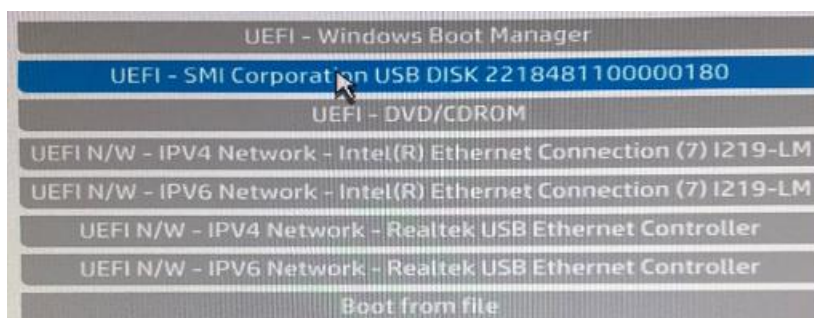
We recommend unplugging any network cable before re-imaging.

If there is an active network connection the re-image process cannot be completed without entering a Microsoft Account. Network connectivity can be established later, also the Microsoft registration can be done later.

- 1 Plug the Recovery USB Flash Drive in the USB port before power-on.
- 2 Press the power button. Immediately press the Escape key repeatedly, about once every second, until the Startup Menu opens.
- 3 Choose **Boot Menu (F9)**



- 4 Choose the USB media, e.g. **UEFI - SMI Corporation USB DISK...**



NOTE:
The actual name may defer from the example in this figure

- 5 The Windows 10 logo will be displayed, followed by a command line window. Enter the command **ApplyImage** there and press ENTER.
- 6 As a next step choose
 - a option 2 if the complete C: drive should be re-imaged as one partition, or
 - b option 1 if a C: and D: partition shall be set up.In doubt choose option 2 (which means no D: partition) because if the user is not aware of the D: drive this could lead to this disk space remaining unused.
- 7 After completing 100% of re-image procedure (see percentage counting) type **exit** to reboot.
- 8 Complete the Windows Setup

NOTE

Windows 10 Pro must be activated as follows:

Once connected to the Internet, the activation may be triggered by Microsoft automatically after some time. Instead waiting for this event, you can actively press [**Windows Key + R**] to open a Run window. Type **slui 4**, and press **<enter>**.

Then choose your country and activate Windows by following the instructions provided.

- 9 ***Restart the Bundle PC to complete the setup script before doing the next steps!!***
(This may take several minutes)

- 10 Double-click the **Adobe** icon on the desktop, accept the license agreement and confirm Adobe Reader as the default PDF application (follow the instructions given > **Apply**)

The PC will now have a unique Device name (*DESKTOP-xxxxxxx*).

Decide if this name should remain or if there are other requirements e.g. from the local IT.

Now proceed with the ChemStation Installation following the instructions provided in the installation manual.

NOTE

For further Microsoft Windows 10 Support, please visit <https://support.microsoft.com>

www.agilent.com

© Agilent Technologies, Inc. 2019

Edition 04/2019
Printed in Germany



M8380-90002

