

Agilent Scan Control V9.1.15 Software Installation Notes

(August 2022)

Scan Control version 9.1.15 is intended for use on Windows 10 64-bit systems. The software was validated with Windows 10 revision 20H2. This note will direct you on how to install Scan Control software version 9.1.15.

Before you begin:

- 1) **Rolling back** the Scan Control version 9.1.15 installation to a previous version (9.1.11 or earlier) of Scan Control requires an Agilent-qualified service engineer.
- 2) If your computer was provided by Agilent it is not necessary to configure the PC network. If you have a non-Agilent supplied PC or you have reimaged the PC, then you may also need to configure the network adapters present in the system to work with the SureScan instrument. See the ***SureScan PC Network Configuration*** instructions included with the Scan Control release.
- 3) If you need to reinstall the current version, select the Repair option when prompted. This will perform a full reinstall.
- 4) Configuration data, such as Scan Regions and Scan Protocols, will be preserved during upgrades and reinstallations. Instrument calibration parameters are stored on the instrument itself and are not affected by reinstallation.
- 5) You must be logged in as an **Administrator** to perform this installation.
- 6) **Managing Windows 10 Updates** - Auto-reboot after a Windows 10 update can interrupt a scanning session. Windows 10 settings can be adjusted to help prevent this. Each revision of Windows 10 provides different controls over updates and auto-reboot. The options available may include deferral of updates (feature and security), setting working hours, delaying updates, or editing the reboot policy. Consult with your IT department to determine the best solution if this is a concern.

To Install Scan Control software, perform the following steps:

- 1) Double click on **Setup.exe** and follow the prompts.
 - Agilent recommends performing installation to the default folder.
 - Setup.exe will first install Agilent Installation Qualification Tool (IQT).
 - It will then install Scan Control software.
 - During the installation, you will be prompted several times by Windows User Access Control to permit installation to proceed. **Please click Yes when prompted.**

Scan Control Installation Notes

- 2) When you first connect to the scanner after installing a Scan Control software update, you will be prompted to install firmware on the instrument.
 - Please follow the instructions presented – you will need to restart the instrument and Scan Control after the firmware is updated.
- 3) After restart, if Scan Control fails to communicate with the scanner, you may need to restore Scan Control Firewall Settings (see Restoration of Firewall Settings below).

Restoration of Firewall Settings

Correct operation of the scanner requires that Scan Control be allowed full access through the Windows Firewall. This rule is set by the installer. If the firewall rules are reset for any reason, you can re-apply the Scan Control rules by executing **SetFirewallRules.bat**, as follows:

- Navigate to the Scan Control executable folder (usually, C:\Program Files (x86)\Agilent\ScanControl)
 - Right-click the file SetFirewallRules.bat
 - Select Run as administrator.
 - Select the allow option.
-