

Release Notes - Scan Control software (August 2022)

This document describes the enhancements and defect fixes related to the release of Scan Control software.

MICROARRAY SCANNER SOFTWARE 9.1.15 (August 2022)

Windows 10 & 11 Rev Compatibility: Scan Control version 9.1.15 is intended for use on Windows 10 & 11 64-bit systems. The software was validated with Windows 10 version 20H2.

Microsoft Windows 10 Automatic Updates: Windows 10 updates can interrupt scanning, see the "Installation Notes" for information on managing Windows 10 updates.

General Software Enhancements:

Fixed Erroneous Laser Over Temperature Errors (Team Track Issues #295950, 272524, 180448, 179756, and 015638). Scan Control software will now detect erroneous laser over temperature errors coming from the Laser Control board for the laser base plate temperature. The erroneous communication error caused false out of range laser temperature errors to be reported to the status system and user. With the new release, if errors are not valid, system will not generate an error message and disrupt operation.

Service Software Enhancements: Improvements and fixes to the "service mode" functionality.

Issue ID: #295890.

MICROARRAY SCANNER SOFTWARE 9.1.13 (January, 2020)

Windows 10 Rev Compatibility: Scan Control version 9.1.13 is intended for use on Windows 10 64-bit systems. The software was validated with Windows 10 version 1903.

Microsoft Windows 10 Automatic Updates: Windows 10 updates can interrupt scanning, see the "Installation Notes" for information on managing Windows 10 updates.

Rolling Back the Installation: Rolling back the Scan Control version 9.1.13 installation to a previous version of Scan Control requires an Agilent qualified service engineer.

General Software Enhancements:

Minimizes cassette movement during barcode reading. Issue ID: 180703

Self-test start requirements identified to user if conditions not met to run self-test. Issue ID: 269931

Error Message and Log File Enhancements

Fixed typos in door closing warning message. Issue ID: 204761

Removed misleading "Activity 'CalibADC' failed" error message from Status Log. Issue ID: 207180

Removed error message when Feature Extraction program path is not set in preference settings. Issue ID: 229285

Improved PMT calibration messages by identifying red or green channel. Issue ID: 229547

Removed erroneous Bar Code Reader Timeout messages from Status Log. Issue ID: 239472

Fixed Scan Control error when Exception Log file is locked by the operating system. Issue ID: 257271

Improved autofocus hold percentage warning statement in Syslog File. Issue ID: 262347

Improved PMT voltage error messages in Syslog File. Issue ID: 265181

Corrected the mismatched laser channel names in the "laser power error" message. Issue ID: 267498

Stopped repeated messages for green laser end of life warning in Status Log. Issue ID: 286550

Scan date/time format changed to 24 hour format for clarity in scan TIFF file. Issue ID: 290200

Service Software Enhancements: Improvements and fixes to the "service mode" functionality.

Issue ID: 159738, 177168, 178252, 178253, 181069, 181688, 183795, 188013, 188178, 190373, 191108, 207646, 213362, 229024, 229026, 229072, 229273, 229282, 229283, 234421, 237288, 239380, 240417, 240424, 250401, 252524, 256594, 258010, 258722, 261315, 262383, 262384, 269928, 269932, 270949, 270950, 271197, 271751, 282002, 283637, 285769, 289594, 291942

MICROARRAY SCANNER SOFTWARE 9.1.11.13 (May 23, 2019)

Windows 10 Rev Compatibility: Scan Control validated to be compatible with Windows 10 rev 1803.

Scan Control User Guide for RUO SureScan G4900DA bundle includes an older version of the User Guide. The latest version of the G4900DA SureScan Scan Control User Guide be accessed via the Agilent.com website. Search: G4900DA User Guide.

MICROARRAY SCANNER SOFTWARE 9.1.11.13

Windows 10 Rev Compatibility: Scan Control updated to provide compatibility with Windows 10 rev 1703. Prevents possible unexpected termination of the software when manipulating application windows. Issue ID 280417

MICROARRAY SCANNER SOFTWARE 9.1.11.7

Error Message Enhancements

Scan Control error messages for fan failures now provide the user with details on which specific fan has failed. Issue ID 229276

Scan Control now provides the user with earlier notice of internal temperature issues that could affect optimal operational performance of the instrument. Issue ID 244710

Scan Control now has enhanced laser power monitoring checks and will report laser power issues both at startup and during normal operation. Issue ID 245248

Operational Safety Enhancement

Scan Control now ensures shutdown of all cassette motion when the cassette door is forced open during the initialization process. Issue ID 244712

Software Regulatory Statement

The "About Agilent Scan Control" box now includes the following statement for the SureScan G4900DA/G2600D model "For Research Use Only. Not for use in diagnostic procedures." Issue ID 253750

Service Software Enhancements

Issue ID 171149, 184853, 189199, 178322, 229022, 229023, 229025, 229275, 229290, 229936, 233401, 236403, 244129, 244140, 244142, 244620, 244688, 244689, 240111, 245436, 247099

MICROARRAY SCANNER SOFTWARE 9.1.9

User Guide Selection

Scan Control now includes a new User Guide selection feature that allows to end user to select the appropriate version of the User Guide to display based upon their scanner model, language, and regional preferences. Issue ID 230537

Service Software Enhancements

Issue ID 020035, 022075, 022111, 233918, 238850

MICROARRAY SCANNER SOFTWARE 9.1.7

Laser end-of-life warning defect fixed

In a few scanners, a "laser end-of-life warning" was displayed in error, due to occasional current variations during laser warm-up. Lasers are now checked after warm-up. Issue ID 195614.

Scan mode added to TIF headers

The TIF header of the image file now indicates if the scan was run in high-sensitivity mode. Issue ID 195966.

Scanning robustness improved for contaminated slides

The auto-focus controls now use additional focus locations to mitigate dirty areas on slides. Issue ID 180200, 195618, 181332.

Additional checks in Self Test

The user Self Test now includes additional checks on laser performance. Issue ID 187593.

Warm-up message

A “Lasers Ready” message was added to the status bar to confirm when lasers have fully warmed up. Issue ID 017362.

Service Software Enhancements

Issue ID 178431, 182974, 183184, 184345, 185202, 186618, 187793, 187795, 189385, 189826, 191318, 198341, 198349, 200378.

MICROARRAY SCANNER SOFTWARE 9.1.5

Service Software Enhancements

Issue ID 181835, 182019, 182085, 182600, 182882.

Installation Qualification Procedural Improvements

Issue ID 182106, 182149, 182752, 182980, 182991.

MICROARRAY SCANNER SOFTWARE 9.1.3

Enhanced sensitivity scanning mode

Scan Control now includes a new, high-sensitivity mode available at all scan resolutions, providing the user with even lower detection limits. Sensitivity is comparable to the double-pass mode but with scan times reduced by 25%. The high-sensitivity mode can be selected in the Resolution pull-down menu of the Scan Protocol Editor. Issue ID 176735.

WARNING: Scan Control 9.1.1 does not recognize the new high-sensitivity mode. After 9.1.3 or a later version is installed, user cannot downgrade back to 9.1.1 unless all high-sensitivity protocols have been deleted or replaced with standard protocols. Issue ID 178924.

Self Test defect fixed

In Scan Control 9.1.1, failure to restart Scan Control after running Self Test can affect the PMT settings and impact the instrument calibration as a result. Restarting the software is no longer necessary in the new 9.1.3 version. Issue ID 176471.

Improved Auto-Focus performance

Auto-Focus performance was improved for non-standard thickness slides to match the precision of standard slide focus. Issue ID 177422.

Slide Queue defects fixed

In rare cases, attempting to add slide slots with blank or custom protocols to the queue to scan could result in unexpected behavior or a software error. The issues were fixed in the new 9.1.3 version. Issue ID 175335, 181313.

In rare sequence of protocol editing selections, PMT gain could be set below its lower limit and trigger an error. Such possibility was eliminated. Issue ID 179753.

WARNING and ERROR statements changed

A “license file not found” WARNING statement was displayed in the log file unless a high-resolution license had been installed, giving the user the impression that something was wrong. Standard resolution scans do not require any license. The warning was modified accordingly for greater clarity. Issue ID 177122.

Error window “folder is not write-accessible” was displayed when user attempted to assign a non-existing scan data folder. The error message was modified to “folder does not exist” for greater accuracy. Issue ID 178980.

If connectivity to the scan data drive is lost during a scan, a new message informs user in which local drive folder the scan was saved. Issue ID 180847.

KNOWN ISSUES

The bottom status bar may not refresh automatically after the PC exits the screensaver mode. Clicking anywhere in the Scan Control window will refresh the status bar.