



Aria Support Tool Release Notes

This document describes the enhancements, defect fixes, and software compatibility issues related to the releases of Aria Support Tool software, which works with both the AriaMx and AriaDx systems.

Aria Support Tool 1.1

Release Date
<ul style="list-style-type: none">• March 2023
Enhancements
<ul style="list-style-type: none">• Analyze block contamination test results and display contaminated wells in a well diagram format and impacted optical module types via existing diagnostic file or from an Aria instrument.• Displays results from pre-requisite diagnostic tests as “pass” or “fail”. Determines if failed block contamination result was due to a failed pre-requisite diagnostic test.• Ability to copy off the diagnostic file from the instrument to a user-designated location. Tool also generates an PDF form of the analysis results and saves to the user-designated location.• Displays revision number in GUI.
Defect Fixes
<ul style="list-style-type: none">• Defect ID: DGGAST-272. Fixed defect where the Motor Calibration verification pass/fail algorithm compares the forward and backward sets separately, against motor calibration specifications.

Aria Support Tool 1.0

Release Date
<ul style="list-style-type: none">• August 2022

Enhancements
<ul style="list-style-type: none"> • User capability to verify motor calibrations as passing or failed via existing filter.txt or from an Aria instrument • Ability to copy off motor calibration files from the instrument to user-designated locations
Defect Fixes
<ul style="list-style-type: none"> • N/A, first release

Known Issues & Workarounds

Known Issues	Workaround
Inaccurate error message shown when PC is not connected to network. The message “Unable to get instrument details” is shown instead of “Unable to connect to instrument” when the Tool is used on a PC that is not connected to any network.	If you receive the above error message “Unable to get instrument details”, please verify your PC’s connection to the network.
Software unable to download the diagnostic file from the instrument if the file name contains a “#” symbol.	Rename file and remove “#” symbol.
Software unable to download the diagnostic file from the instrument if the file name contains 225 characters.	Reduce number of characters in file name.