Aria Software Revision History and Release Notes

Revision	Release Date	Enhancements, Fixes, Known Defects, and Workarounds
2.0	Release Date September 2022	 Enhancements, Fixes, Known Defects, and Workarounds Enhancement: Ability to verify that the data incoming from an Aria instrument is coming from the matching source. AriaDx Firmware [AriaDx_Firmware_3_1_2202_0902_FPR] Enhancement: Include instrument identifier information in data transmitted to the software. AriaMx Firmware [AriaMx_Firmware_3_1_2202_0901_FPR] Enhancement: Include instrument identifier information in data transmitted to the software. AriaMx Firmware [AriaMx_Firmware_3_1_2202_0901_FPR] Enhancement: Include instrument identifier information in data transmitted to the software. Compatibility Notice: Please note that this AriaDx firmware v2.0 (3.1.2202.0902) and AriaMx firmware v2.0 (3.1.2202.0901) are compatible with software v2.0 (2.0.2204.0441) and above ONLY. There is a risk of the data mismatch issue occurring when using a software release version earlier than version 2.0 to monitor an experiment run on firmware version 2.0. Known Defects & Workarounds Note: These defects will not impact data collection or storage Defect #1 – Users who 1) run the Electronic Tracking (ET) software, 2) run a User Defined Experiment template, and 3) monitor the run could have their SW crash at the end of the run. All 3 conditions must apply and there is no threat to losing data. Workaround #1 – Relaunch the Aria SW to continue data analysis. Defect #2 – Users who 1) "Generate Report" from the Aria SW and 2) have their Windows Display Settings set to 125% or above will have the scaling of the resulting .pdf report reduced.
		 Workaround #2 – Close the Aria SW application and change the scaling of the Windows Display Settings to 100%. Then re-open the Aria SW and re-generate the report.
1.8	March 2021	 Enhancement: Ability to automatically hide the low signal "noise" in the log plots view Enhancement: Support for the import/export of LIMS files in csv format Fixed: Analysis Term Setting: Amplification Plot Smoothing GUI not updated when trigger run from post file Fixed: Analysis Term Settings: Threshold fluorescence not updated when reopening an experiment file Fixed: Analysis Term Settings: Cycle range not applied when importing a thermal profile, then switching to a new tab after first data collection Fixed: Analysis Term Settings: Baseline Correction is not applied for a new run when current active tab is not at monitoring run tab AriaDx Firmware: [AriaDx_Firmware_3_1_2012_2202_FPR]

		 Enhancement: Allow for instrument warmup Fixed Consists are executed an execution discusses to state
		 Fixed: Genesis.exe error when performing diagnostic tests
		 Fixed: Instrument warm up message not appearing after
		instrument health check
		AriaMx: [AriaMx_Firmware_3_1_2012_2201_FPR]
		• Enhancement: Allow for instrument warmup
		• Fixed: Genesis.exe error when performing diagnostic tests
		• Fixed: Instrument warm up message not appearing after
		instrument health check
		Compatibility notice: Please note that this AriaDx firmware
		(3_1_2012_2202) and AriaMx firmware (3_1_2012_2201) are compatible
		with software release versions 1.8 and 1.7.1 ONLY. Experiment files with
		new thermal profile settings (multiple data collection within an
		Amplification segment or Touch down PCR settings) created by this
		firmware version is not supported by software release versions prior to
		1.7.1.
1.7.1	November 2019	Corrected defect discovered in version 1.2 and later which could prevent
		Cq values being included in the Results Table.
1.7	April 2019	Thermal profiles allow multiple data collection points on different
		plateaus in the amplification segment
		• Time and temperature increment settings added to amplification segment
		in thermal profile setup, which allows for Touch Down PCR
		In Relative Quantity Chart, now we will be able to show that selected
		wells have multiple targets
		Added ability to assign amplification plot colors by dye or target
		Plate Setup update: Now, sample names remain unchanged when the
		target names are changed
		Adaptive Baseline correction: Consistent with prior Aria software release
		versions 1.3, 1,4, and 1.51, the Adaptive Baseline correction is now the
		default setting in Aria software release ver. 1.7.
1.6	November 2017	Import well names from Excel files
		Import thermal profiles from existing files
		Display Cq values on the analysis screen
		Keep row sorting when exporting the Result Table to Excel file
		Export chart data in horizontal format
		Save graph settings
		Show data in Standard Curve plot when Standard well type is filtered
		Correct baseline setting in an analysis template
1.5.1	July 2017	Corrected defects discovered in version 1.5
1.5	June 2017	Configure more segments, plateaus and amplification cycles during the
		thermal profile setup
		Save default analysis settings as a template
		Generate a report after each validation plate run
		 Perform multiple column sorting on tabular data

		Export data based on column or row option
1.4.1	May 2017	Corrected several defects discovered in version 1.4
1.4	November 2016	Added LIMS support
1.3	August 2016	 Introduced compatibility with Windows 10, 64bit (32 bit not tested) Fixed a defect that disabled capture of changes to the report files Mx/Dx mode implementation
1.2	April 2016	Added the requirement for License Activation to use HRM analysis
1.1	November 2015	 Implemented crosstalk reduction between spectrally similar optical cartridges to allow for release of Atto425 cartridge, resolving cartridge overlap with FAM. Fixed 2 defects related to the Electronic Tracking (ET) software
1.0	September 2014	Aria first release