

## Aria Software Revision History and Release Notes

Revision	Release Date	Enhancements, Fixes, Known Defects, and Workarounds
2.1.1	September 2024	<ul style="list-style-type: none"> <li>• Enhancement: Support for Windows 11 Pro (64 bit).</li> <li>• Enhancement: Support for Microsoft Visual C++ 2022 Redistributable package (x86).</li> <li>• Enhancement: Support for Microsoft .NET Framework 4.8 installer.</li> <li>• Enhancement: Corrected wording from “CFR” to “ET”.</li> <li>• Fixed: Software no longer crashes while browsing the file explorer when the instrument has a user account name like the instrument name.</li> <li>• Fixed: User can now copy and delete experiment files from instrument to the software in ET mode via the instrument file explorer.</li> <li>• Fixed: The display layout now considers the maximum number of characters supported for the instrument name and username so that a long login username is not truncated on the Instrument Information dialog box.</li> <li>• Fixed: Addressed security vulnerability that’s present in the scenario when the user is re-monitoring an experiment that’s currently running.</li> <li>• Fixed: Updated Technical Support’s contact email</li> <li>• Compatibility Notice: In all cases, it’s always best practice when updating software, to also update firmware <b>if</b> a new revision is available and compatible. Please note that there is no new release of AriaDx firmware or AriaMx firmware with SW v2.1.1. For full performance, run Aria SW v2.1.1 with Aria firmware v2.1 (Dx: 3.1.2306.0601 and Mx: 3.1.2306.0602). Note that this AriaDx firmware v2.1 (3.1.2306.0601) and AriaMx firmware v2.1 (3.1.2306.0602) are compatible with software v2.0 (2.0.2204.0441) and above <b>ONLY</b>. <ul style="list-style-type: none"> <li>○ There is a risk that well signal may be shifted to an adjacent well if the experiment is run after a motor calibration is performed on a firmware release version earlier than v2.1.</li> <li>○ 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 2.1.</li> </ul> </li> </ul> <p><b>Known Defects &amp; Workarounds</b></p> <ul style="list-style-type: none"> <li>• Microsoft Defect #1 – On Win11 PC, when installing Aria SW in Electronic Tracking mode, the installer displays an error message of “Could not connect to Microsoft SQL Server” or “Windows could not start the SQL Server on Local Computer”.</li> <li>• Microsoft Workaround #1 – Refer to troubleshooting guidance from Microsoft, available at <a href="#">Troubleshooting operating system disk sector size greater than 4 KB - SQL Server   Microsoft Learn</a>. You may need to work</li> </ul>

		<p>with your organization's Information Technology group and/or Microsoft Technical Support to resolve the issue.</p> <ul style="list-style-type: none"> <li>• Defect #2: In ET mode, the user is unable to copy a template file off the instrument.</li> <li>• Workaround #2: Copy the template file from the instrument using a USB drive.</li> </ul>
2.1	November 2023	<ul style="list-style-type: none"> <li>• Enhancement: SQL wrapper installer now installs SQL Server 2019 Express Edition for ET SW.</li> <li>• Fixed: SQL wrapper installer now prevents users from installing over an existing SQL server instance.</li> <li>• Fixed: Rephrased the error message when the user clicks "No" to cancel the installation from the "User Access Control" window.</li> <li>• Fixed: Broken "Cancel" button is removed from the installation progress bar.</li> <li>• Fixed: The software now prompts the user to close some experiments when the memory consumption exceeds 1.6GB.</li> <li>• Fixed: On the "My Template" page, the "Database Templates" table now displays the Experiment Type for User-Defined experiments correctly.</li> <li>• Fixed: In ET mode, "Save As" now works correctly when the user right-clicks other tabs while a post-run file is opened on the first tab.</li> <li>• Fixed: On the "Generate Report" page, in the "Report Configuration" panel, the Items list has a scrollbar to prevent the "Generate Report" button from disappearing.</li> <li>• Fixed: In ET Mode, the "Generate Report" button displays correctly when the experiment contains 5 plots.</li> <li>• AriaDx Firmware [AriaDx_Firmware_3_1_2306_0601_FPR] <ul style="list-style-type: none"> <li>○ Enhancement: Implemented verification of motor calibration values against specifications.</li> <li>○ Fixed: The firmware now allows only FAM optical module to be used for motor calibrations.</li> <li>○ Fixed: Removed visual glitch when dismissing a motor calibration error message.</li> <li>○ Fixed: Removed "AriaMx" from the Help File for AriaDx mode.</li> <li>○ Fixed: Message box title now matches the instrument mode.</li> </ul> </li> <li>• AriaMx Firmware [AriaMx_Firmware_3_1_2306_0602_FPR] <ul style="list-style-type: none"> <li>○ Enhancement: Implemented verification of motor calibration values against specifications.</li> <li>○ Fixed: The firmware now allows only FAM optical module to be used for motor calibrations.</li> <li>○ Fixed: Removed visual glitch when dismissing a motor calibration error message.</li> </ul> </li> </ul>

		<ul style="list-style-type: none"> <li>○ Fixed: Message box title now matches the instrument mode.</li> <li>● Compatibility Notice: In all cases, it's always best practice when updating software, to also update firmware if a new revision is available and compatible. Please note that this AriaDx firmware v2.1 (3.1.2306.0601) and AriaMx firmware v2.1 (3.1.2306.0602) are compatible with software v2.0 (2.0.2204.0441) and above ONLY.       <ul style="list-style-type: none"> <li>○ There is a risk that well signal may be shifted to an adjacent well if the experiment is run after a motor calibration is performed on a firmware release version, earlier than v2.1.</li> <li>○ 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 2.1.</li> </ul> </li> </ul> <p><b>Known Defects &amp; Workarounds</b></p> <ul style="list-style-type: none"> <li>● NA</li> </ul>
2.0	September 2022	<ul style="list-style-type: none"> <li>● Enhancement: Ability to verify that the data incoming from an Aria instrument is coming from the matching source.</li> <li>● AriaDx Firmware [AriaDx_Firmware_3_1_2202_0902_FPR]       <ul style="list-style-type: none"> <li>○ Enhancement: Include instrument identifier information in data transmitted to the software.</li> </ul> </li> <li>● AriaMx Firmware [AriaMx_Firmware_3_1_2202_0901_FPR]       <ul style="list-style-type: none"> <li>○ Enhancement: Include instrument identifier information in data transmitted to the software.</li> </ul> </li> <li>● 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.</li> </ul> <p><b>Known Defects &amp; Workarounds</b></p> <p><i>Note: These defects will not impact data collection or storage</i></p> <ul style="list-style-type: none"> <li>● 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.</li> <li>● Workaround #1 – Relaunch the Aria SW to continue data analysis.</li> <li>● 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.</li> <li>● 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.</li> </ul>

1.8	March 2021	<ul style="list-style-type: none"> <li>• Enhancement: Ability to automatically hide the low signal "noise" in the log plots view</li> <li>• Enhancement: Support for the import/export of LIMS files in csv format</li> <li>• Fixed: Analysis Term Setting: Amplification Plot Smoothing GUI not updated when trigger run from post file</li> <li>• Fixed: Analysis Term Settings: Threshold fluorescence not updated when reopening an experiment file</li> <li>• Fixed: Analysis Term Settings: Cycle range not applied when importing a thermal profile, then switching to a new tab after first data collection</li> <li>• Fixed: Analysis Term Settings: Baseline Correction is not applied for a new run when current active tab is not at monitoring run tab</li> <li>• AriaDx Firmware: [AriaDx_Firmware_3_1_2012_2202_FPR]             <ul style="list-style-type: none"> <li>○ Enhancement: Allow for instrument warmup</li> <li>○ Fixed: Genesis.exe error when performing diagnostic tests</li> <li>○ Fixed: Instrument warm up message not appearing after instrument health check</li> </ul> </li> <li>• AriaMx: [AriaMx_Firmware_3_1_2012_2201_FPR]             <ul style="list-style-type: none"> <li>○ Enhancement: Allow for instrument warmup</li> <li>○ Fixed: Genesis.exe error when performing diagnostic tests</li> <li>○ Fixed: Instrument warm up message not appearing after instrument health check</li> </ul> </li> <li>• 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.</li> </ul>
1.7.1	November 2019	<ul style="list-style-type: none"> <li>• Corrected defect discovered in version 1.2 and later which could prevent Cq values being included in the Results Table.</li> </ul>
1.7	April 2019	<ul style="list-style-type: none"> <li>• Thermal profiles allow multiple data collection points on different plateaus in the amplification segment</li> <li>• Time and temperature increment settings added to amplification segment in thermal profile setup, which allows for Touch Down PCR</li> <li>• In Relative Quantity Chart, now we will be able to show that selected wells have multiple targets</li> <li>• Added ability to assign amplification plot colors by dye or target</li> <li>• Plate Setup update: Now, sample names remain unchanged when the target names are changed</li> <li>• 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.</li> </ul>
1.6	November 2017	<ul style="list-style-type: none"> <li>• Import well names from Excel files</li> </ul>

		<ul style="list-style-type: none"> <li>• Import thermal profiles from existing files</li> <li>• Display Cq values on the analysis screen</li> <li>• Keep row sorting when exporting the Result Table to Excel file</li> <li>• Export chart data in horizontal format</li> <li>• Save graph settings</li> <li>• Show data in Standard Curve plot when Standard well type is filtered</li> <li>• Correct baseline setting in an analysis template</li> </ul>
1.5.1	July 2017	<ul style="list-style-type: none"> <li>• Corrected defects discovered in version 1.5</li> </ul>
1.5	June 2017	<ul style="list-style-type: none"> <li>• Configure more segments, plateaus and amplification cycles during the thermal profile setup</li> <li>• Save default analysis settings as a template</li> <li>• Generate a report after each validation plate run</li> <li>• Perform multiple column sorting on tabular data</li> <li>• Export data based on column or row option</li> </ul>
1.4.1	May 2017	<ul style="list-style-type: none"> <li>• Corrected several defects discovered in version 1.4</li> </ul>
1.4	November 2016	<ul style="list-style-type: none"> <li>• Added LIMS support</li> </ul>
1.3	August 2016	<ul style="list-style-type: none"> <li>• Introduced compatibility with Windows 10, 64bit (32 bit not tested)</li> <li>• Fixed a defect that disabled capture of changes to the report files</li> <li>• Mx/Dx mode implementation</li> </ul>
1.2	April 2016	<ul style="list-style-type: none"> <li>• Added the requirement for License Activation to use HRM analysis</li> </ul>
1.1	November 2015	<ul style="list-style-type: none"> <li>• Implemented crosstalk reduction between spectrally similar optical cartridges to allow for release of Atto425 cartridge, resolving cartridge overlap with FAM.</li> <li>• Fixed 2 defects related to the Electronic Tracking (ET) software</li> </ul>
1.0	September 2014	<ul style="list-style-type: none"> <li>• Aria first release</li> </ul>