



Addendum to AriaMx Manual

Tips for updating your firmware and PC GUI software

This communication will highlight run a software update on the AriaMx Real-Time PCR System.

(1) Go to the website:

www.agilent.com/genomics/ariamxsoftware

This is a vanity website, so it actually redirects. You will see this screen:

The screenshot shows a web browser window with the following elements:

- Address bar: <http://www.genomics.agilent.com/article.jsp?pageId=56000048&CID=G011530>
- Navigation menu: File, Edit, View, Favorites, Tools, Help
- Search bar: Find: malaysia
- Page header: PRODUCTS & SERVICES SOLUTIONS
- Main content area:
 - AriaMx Software Download**
 - AriaMx Software (instrument control and data analysis)
 - Download AriaMx PC Software Update: Revision v1.0.1408.2642
 - AriaMx Instrument Firmware
 - No updates currently available

(2) For updating PC software, download to a your PC.

The description of the bug and process is below:

PRODUCTS & SERVICES SOLUTIONS TECHNICAL SUPPORT

AriaMx PC Software 1.0 Update

Note: The AriaMx PC Software is the user interface software for PC that interfaces with the instrument. This software is usable only with the Agilent AriaMx (G8830A). It will not run on Linux systems.

AriaMx PC Software Update: Revision 1.0.1408.2642

Revision 1.0.1408.2642 was released in September 2014 to correct a defect where selecting the ctrl key in the Raw Data Plots page could cause the application to close unexpectedly. This update resolves the following issues which could cause the application to close:

- Opening a corrupted experiment after successfully opening an experiment and project on same tab
- Attempting to run an experiment using a previously saved pre-run file, after successfully opening an experiment and project on same tab

Download and Install Instructions

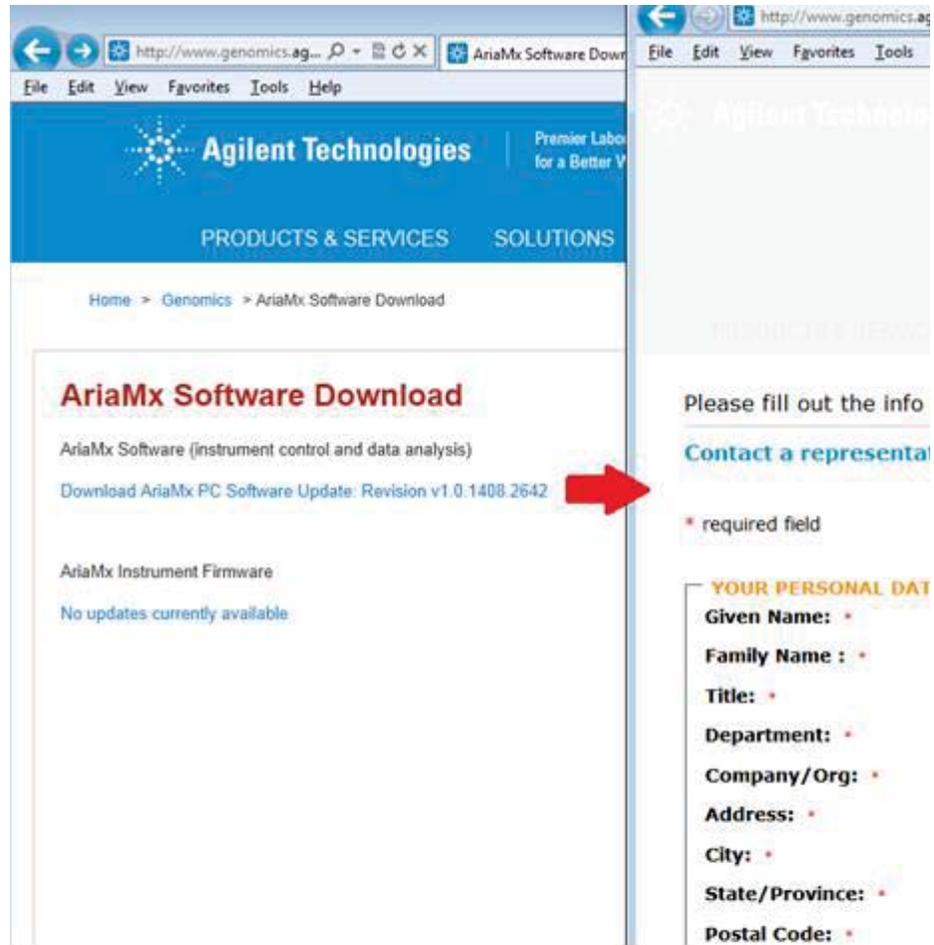
1. [Download](#) the AriaMx PC Software 1.0 Update file (ZIP, 25Mb) to your PC.
2. Unzip the downloaded file to create folder "AriaMx PC Software 1.0 Installer".
3. Manually uninstall any AriaMx PC software already installed on the PC. This can be done from the Control Panel found in the pop-up panel that appears when the Start button (lower left) is clicked. Choose *Control Panel > Programs > Programs and Features* option to display the Control Panel. Find "Agilent AriaMx 1.0" in the list, right-click on it and then click **Uninstall**.
4. Follow the steps described in the [Installation of the AriaMx Software](#) section of the [Setup and User Guide](#) to install the PC software (using the folder created in step 2 above rather than the default folder).

Note: If installing the ET (Electronic Tracking) software, you must first have the SQL Server Instance installed as described in the [Setup and User Guide](#). If not yet installed, the SQL Server Instance software CD delivered with the instrument.

If you have trouble installing your software update, you can [contact technical support](#)

Please note you will need to enter personal information after clicking on the "Download AriaMx PC Software Update" link in order to get to the actual download page.

During the actual software installation you will need to accept a license agreement.



(3) For updating firmware

- a. Wait until Jan 2015
- b. Go to the same website as above
- c. Click the download link under firmware
- d. It is likely you will have to download this to a USB stick
- e. Log in as an administrator to the AriaMx instrumentation
- f. Install software from USB to AriaMx.
- g. Should take about 8 to 10 mins.

WHERE TO FIND RESOURCES

Technical support:

Genomics Division- Bioreagents team
800-227-9770-3-4-3 tel.
858 777 5387 fax

Collaborations: Please contact Mary Napier, Senior Product Manager Genomics and Molecular Diagnostics, or Caroline Tsou if you have any questions regarding this product.

Software updates: www.agilent.com/genomics/ariamxsoftware



For Research Use Only. Not for use in diagnostic procedures.