

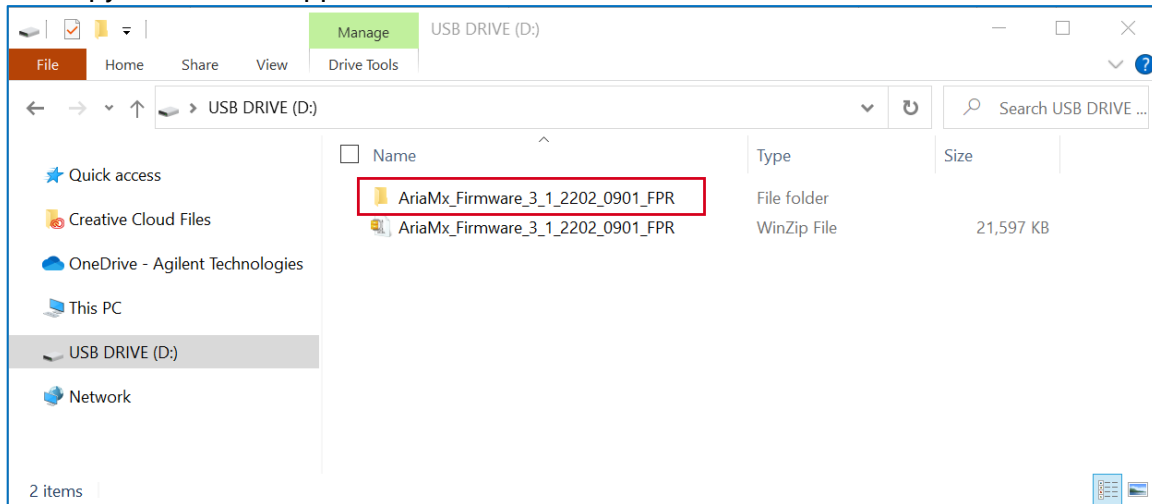
AriaMx/Dx How to upgrade or load instrument firmware

Prepare Firmware (FW)

1. Download the latest FW revision from Agilent onto PC.

Note: AriaMx FW must only be installed on AriaMx Instruments and AriaDx FW must only be installed on AriaDx instruments. The latest Aria FW revisions may be found on the Aria Software and Firmware download pages. FW will download as a .zip file.

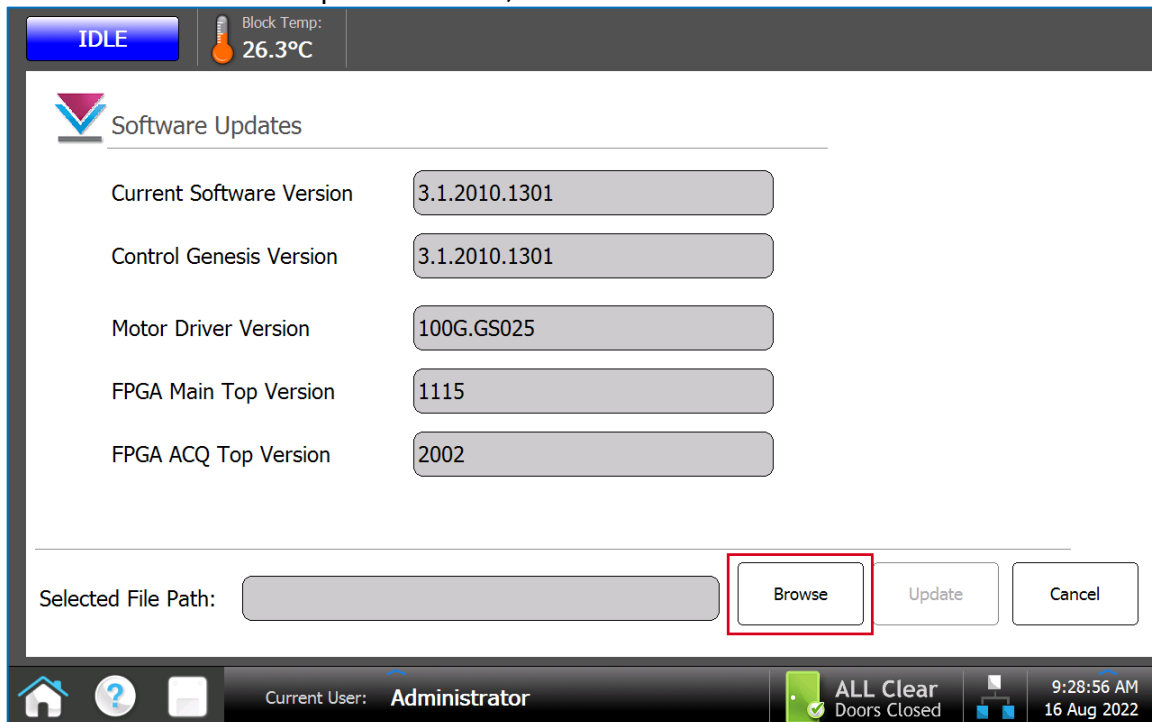
2. Copy and save unzipped FW folder onto a USB Drive



3. Insert USB Drive with FW into the front or rear USB port of the Aria Instrument

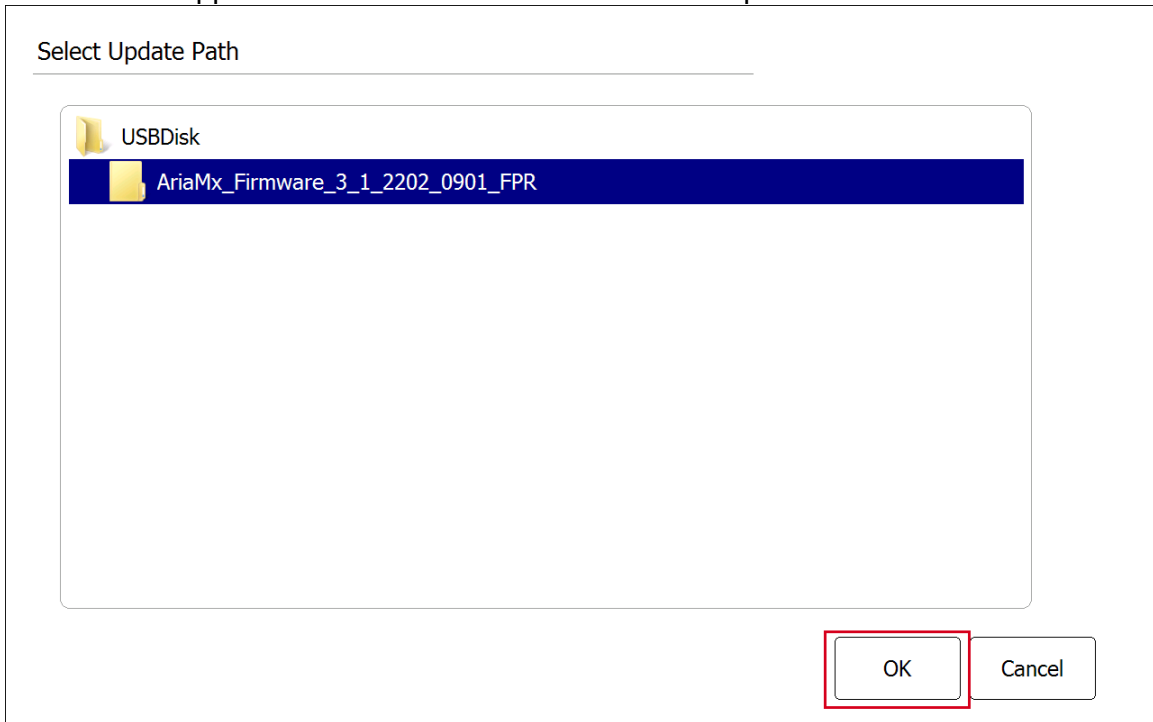
How to upgrade/load instrument firmware

1. From the Instrument Home Screen, log in as **Administrator**
2. From the instrument **Home Screen** >> **Settings** >> **Software Updates**
3. From the Software Updates Screen, Select **"Browse"**

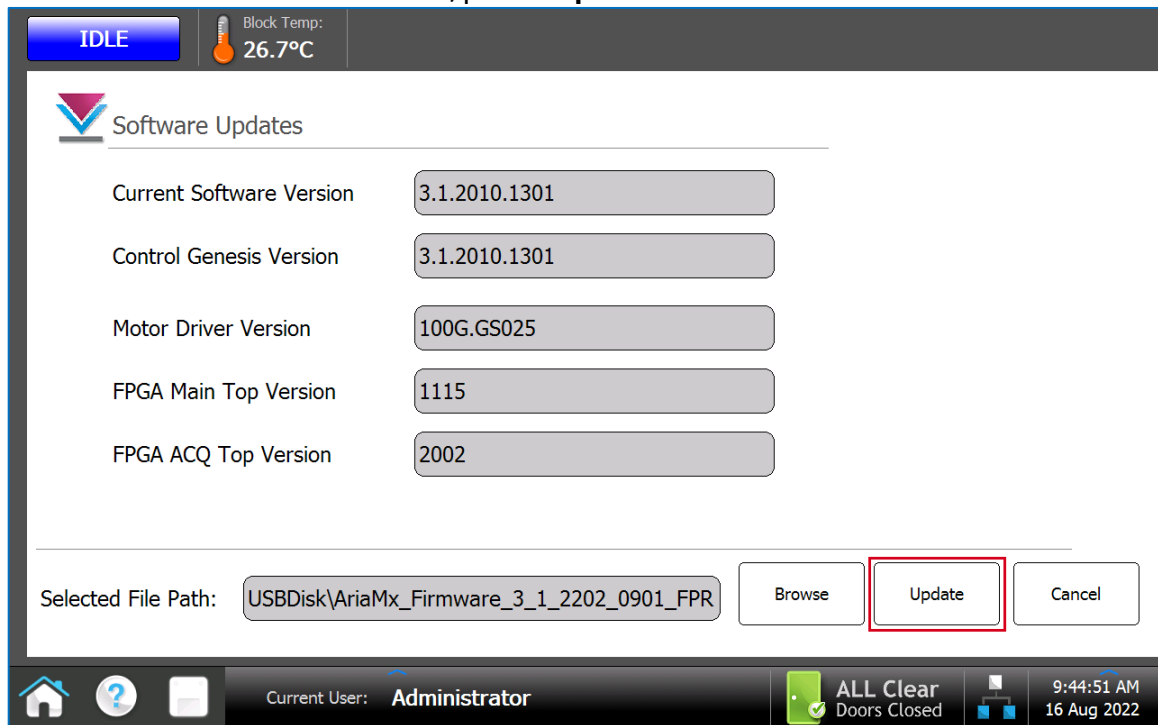


4. Double-click on **"USB Disk"** folder

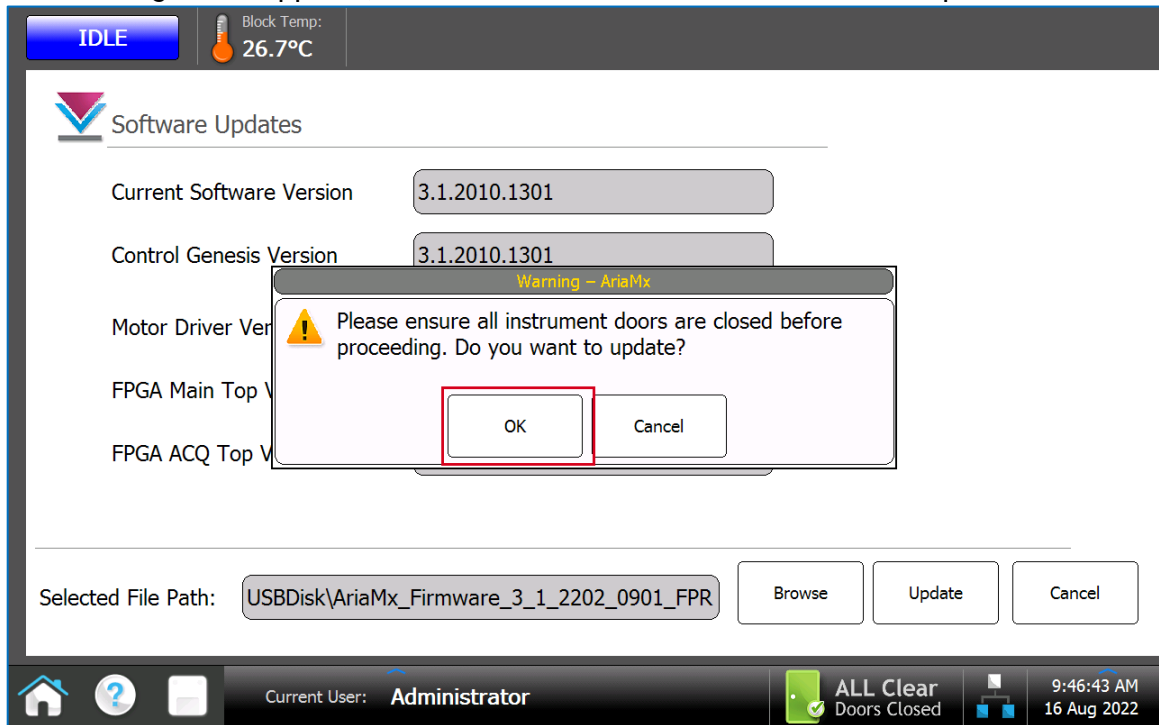
5. Select the unzipped Firmware folder on the USBDisk and press "OK"



6. With the correct File Path selected, press "Update"



7. When Dialogue box appears, ensure all instrument doors are closed and press "OK"



8. After Instrument firmware has been uploaded, Aria instrument will automatically reboot.
Note: It is critical that the Aria instrument is rebooted for instrument firmware upgrade to successfully complete. If Aria Instrument does not automatically reboot after FW upgrade, manually reboot Aria Instrument.

End