

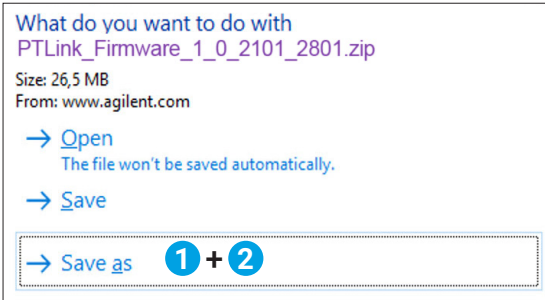
PT200 Link

Steps of Updating Firmware

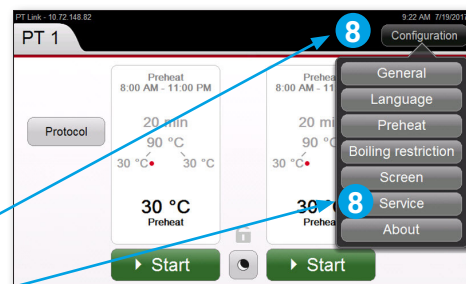
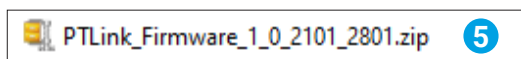
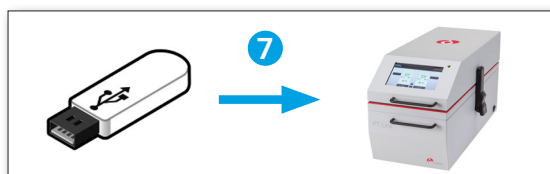
Please check your PT Link(s) (PT20027 and/or PT20040) to verify that it has the most recent software update, version: PTLINK_Firmware_1_0_2101_2801.

This information can be viewed from the "About screen". To get to the "About" screen, select the 'About' option from the "Configuration" menu. If additional information is needed to check the version number please reference "About PT Link" in the user guide section 7, also called "Configuration Options".

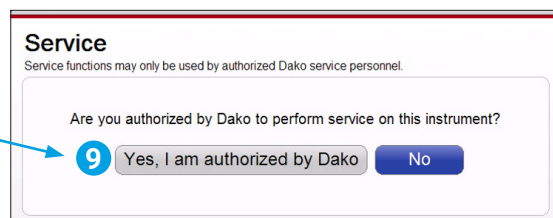
1. Download the PT200 Firmware update file from:
http://www.agilent.com/cs/library/software/public/PTLink_Firmware_1_0_2101_2801.zip
2. Click "save as" to save the zip-file in a folder on your computer
3. Click "open" to unzip the zip-file
4. Make sure the zip-file contains at least these files:
5. **Copy** the folder containing the files (as shown in the picture to the right) to a USB drive
Note: USB drive should have at least 40 MB of free storage, and needs to be **compatible with** the USB port on the PT Link (USB Type A)
6. Make sure the folder on the USB contains at least these files:
7. Remove the USB from your computer by safely ejecting it and plug it in to the USB port on the PT200 instrument
Note: If you need to move the PT200 instrument to access the USB drive located on the back, please make sure to remove the PT200 tanks to avoid spilling the Target Retrieval Solution
8. Click **Configuration** on PT200, and then click **Service**



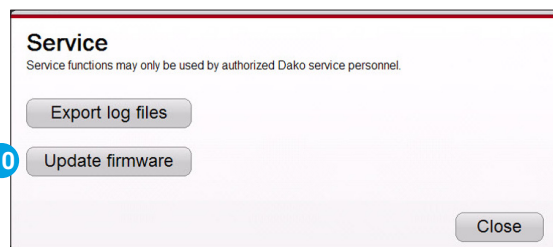
| Name | Type | Size |
|--------------------|--------------|-----------|
| BootImage_B0.nb0 | NB0 File | 768 KB |
| FontsInstaller.CAB | WinZip File | 14,337 KB |
| index.xml | XML Document | 1 KB |
| ptlink_top.bin | BIN File | 1,450 KB |
| xip.bin | BIN File | 33,893 KB |



9. Click **Yes, I am authorized**

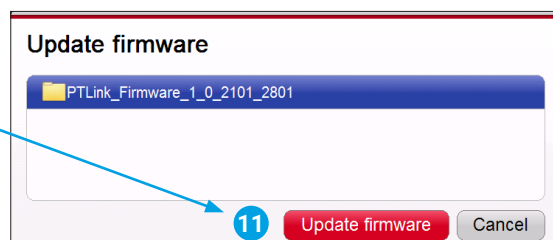


10. Click **Update firmware**

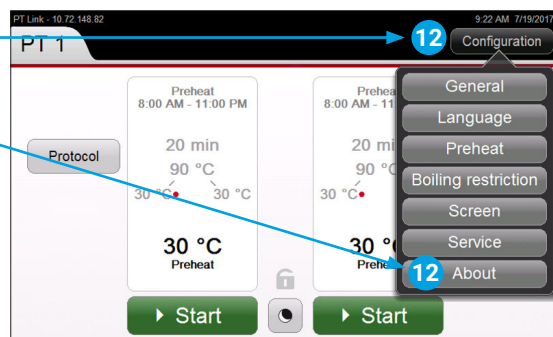


11. Select the firmware to be updated, and then click **Update firmware**

Wait until the update is completed successfully and the instrument is **rebooted automatically**



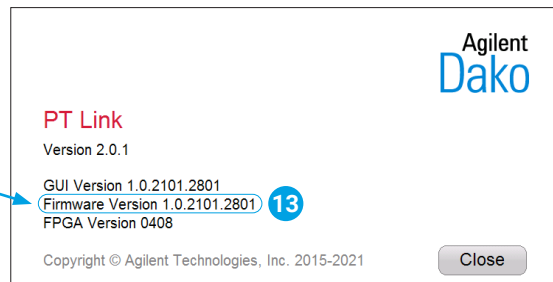
12. Click **Configuration**, and then click **About**



13. Check if the installed **instrument firmware version matches the downloaded version**. In this example, the firmware version shall be 1.0.2101.2801.

14. Firmware update completed

15. Remove USB from the USB drive on the PT200



www.agilent.com

This information is subject to change without notice.