
Installation Notes: FFA 120 v2 Assay file (Rev A)

Installation procedure on LabChip® GX/GXII Touch™ instruments:

1. Check the software version that is currently installed on your LabChip® GX Touch™ or LabChip® GXII Touch™ instrument. If it is lower than version 1.10, it is necessary to be updated to version 1.10 or higher.
Please find the latest LabChip® GX Touch™ Software (Not GX Reviewer) from the following link: <https://www.revvity.com/software-downloads/labchip> and follow the instructions provided for software version update.
2. Close the LabChip® GX Touch™ Software.
3. Download the assay file “FFA120 v2.asyx” from above software downloads website.
4. Copy the assay file “FFA120 v2.asyx” into the assay folder “C:\...\LabChip GX Touch\Assay)” on the computer of LabChip® GX Touch™ or LabChip® GXII Touch™ instrument.
5. Reopen the LabChip® GX Touch™ Software.
6. Load the correct LabChip (DNA 5K/RNA/CZE HT LabChip), and then select **HT small RNA** as the chip type from the pop-up window.
7. Confirm the “FFA120 v2” assay is available in the assay drop-down list and select it.

Installation procedure on PC which is not connected to LabChip® GX/GXII Touch™ instruments for data review:

1. Use LabChip® GX Reviewer Software to review data created by “FFA 120 v2.asyx”. “FFA 120 v2.asyx” installation is ONLY for instruments’ operation; it is not required for data reviewing or data analysis with LabChip® GX Reviewer Software.
2. If the LabChip® GX Reviewer Software version is lower than 5.8, please update it to version 5.8 or above. Please find the latest LabChip® GX Reviewer Software from the following link: <https://www.revvity.com/software-downloads/labchip> and follow the instructions provided for software version update.