

Program Update Instructions for Amaxa® Nucleofector® II Device

To view the current software version that comprises the operating programs of your Amaxa® Nucleofector® II Device, turn on your device first (orange button). Press “menu” repeatedly until menu item “Software versions” (for details please refer to the Nucleofector® Manual provided with your device). If the third number displayed is S3-8, your device does not require an update. If it is S3-7, or lower an update is required. You can easily update the programs by using the Amaxa® Chip Card and following the instructions below.

Instructions



1. Turn the device on (orange button).
2. In order to upgrade* the program list in your Nucleofector® II Device, insert the chip card (chip forward, facing down) into the chip card reader slot (see picture).

Do not remove the chip card from the slot before the entire process is finished.

3. After detecting the chip card, the display shows “Card found” and “Read chipcard?”. Press the “X” button to confirm.
4. A running bar indicates the reading of the chip card. After doing so “Success” and “Pulstable 2 stored” is indicated. Press the “X” button.
5. Now the display shows “Progr. found” and “Load programs?”. Press the “X” button again.
6. After a running bar has indicated the loading of the programs the display shows “OK” and “Remove card”.
7. Remove card and press the “X” button twice.
8. Press “menu” repeatedly until menu item “Software versions”. Now the third number displayed is S3-8. The device is updated.

* The stored custom programs will be neither deleted nor overwritten by upgrading novel non-custom optimized programs.