

## Program Update Instructions for Amaxa® Nucleofector® I Device

---

To find out whether it is necessary to update the programs on your Nucleofector® I Device, you can check the software version of your device by simply switching it on. If the second number presented on the display is V2.5, your device does not require an update. If the number is **not** V2.5, you can easily update the programs by using the Amaxa® Chip Card and following the instructions below.

### Instructions

---



1. Turn on the Nucleofector® Device.
2. Insert the chip card (chip forward and downward) into the small slot below the control panel within the grip of the Nucleofector® I Device (see picture).
3. To load the new programs press the “X” button.

---

A successful update is indicated as “C OK” shown on the display.

---

4. Remove the chip card from the Nucleofector® Device
5. Confirm by pressing the “X” button.

---

The display will first show either “R2.0”, “R2.1”, “R2.2”, “R2.3” or “R2.4” (depending on the age of the Nucleofector® Device) and shortly afterwards V2.5.

---

---

The Nucleofector® Device is now updated and ready for use.

---