

AMPLIFICATION

Checking for Firmware and Software Updates for the MiniOpticon™ Real-Time PCR Detection System — A Quick Guide

Bio-Rad Laboratories, Inc. periodically offers updates for the MiniOpticon system firmware and CFX Manager™ software. We strongly recommend that you regularly check your system and install any available updates. Installing updates ensures optimal performance from your instrument and provides additional features as they become available.

To confirm that your system is up to date, check the version of firmware installed on the MiniOpticon system and the version of CFX Manager software installed on your computer. Compare these to the latest versions posted on the Bio-Rad website at www.bio-rad.com/pcr_updates/.

Checking the Firmware Version Installed on the MiniOpticon System

1. Turn on the instrument; wait until the MiniOpticon system is fully booted and shows the main menu screen.
2. In the front panel of the MiniOpticon system, use the arrow keys to navigate to **TOOLS** in the main menu (Figure 1), then press **Enter**.
3. In the Tools menu, navigate to the **VERSION** menu item (Figure 2), then press **Enter**.
4. Compare the installed firmware version displayed on the instrument (Figure 3) with the most current firmware version available on the website.
5. If the version installed on your instrument does not match the current version listed on the website, download and install the update.



Fig. 1. Main menu.

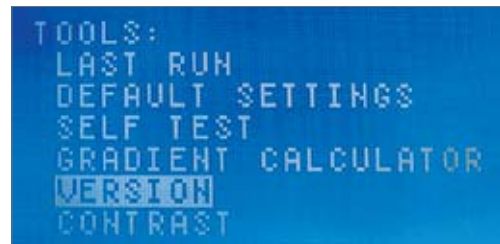


Fig. 2. Tools menu.

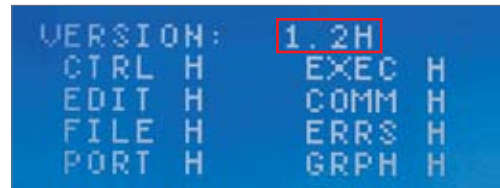


Fig. 3. Installed firmware version displayed on the MiniOpticon system.

Checking the CFX Manager Software Version Installed on Your Computer

1. Launch CFX Manager software.
2. Select **Help > About** from the main software window (Figure 4).
3. The About Bio-Rad CFX Manager window displays the installed software version (Figure 5).
4. If the version installed on your computer does not match the current version listed on the website, download and install the update.

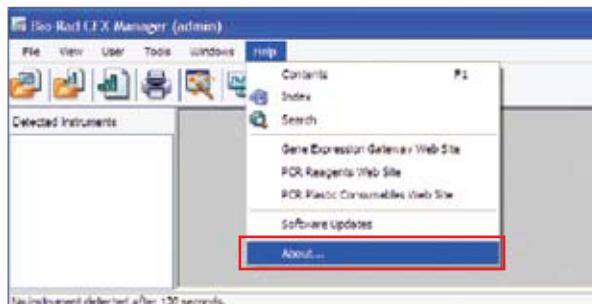


Fig. 4. Navigating in the main window of CFX Manager software.



Fig. 5. Installed software version displayed in the About Bio-Rad CFX Manager window.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applera's United States Patent 6,814,934 B1 for use in research, human in vitro diagnostics, and all other fields except veterinary diagnostics.

Bio-Rad's real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Nos. 6,767,512 and 7,074,367.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400
Canada 905 364 3435 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300 Israel 03 963 6050
Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000