

AMPLIFICATION

Checking Firmware and Software Updates for the C1000 Touch™ Thermal Cycler and the CFX96 Touch™ and CFX384 Touch™ Real-Time PCR Detection Systems — A Quick Guide

Bio-Rad Laboratories, Inc. periodically offers updates for C1000 Touch thermal cycler, CFX96 Touch system, and CFX384 Touch 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 C1000 Touch thermal cycler and the CFX96 Touch or CFX384 Touch 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 C1000 Touch Thermal Cycler and the CFX96 Touch or CFX384 Touch System

1. Turn on the instrument; wait until the system is fully booted and shows the main menu screen.
2. On the main menu screen of the C1000 Touch thermal cycler, touch the **Tools** button (Figure 1).
3. Touch **About** in the User menu (Figure 2).
4. Compare the installed firmware version displayed on the instrument (Figure 3) with the most current firmware version available on the website.

Note: The 8051 and DSP firmware are only used for real-time PCR. If you are using a C1000 Touch thermal cycler for standard PCR, these firmware versions will not be visible.

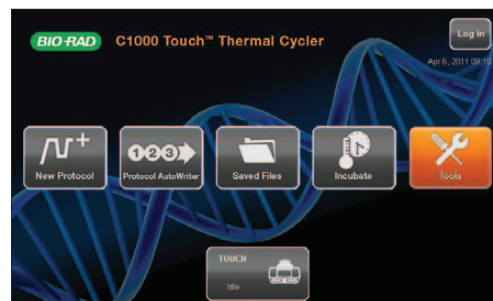


Fig. 1. Main menu of the C1000 Touch thermal cycler.

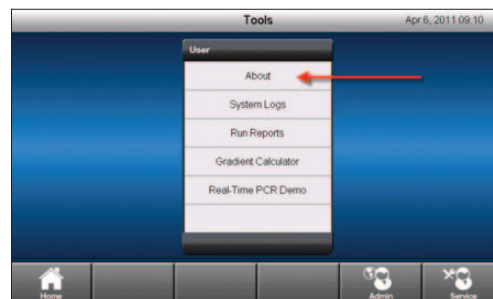


Fig. 2. User menu.



Fig. 3. Installed firmware versions displayed on the C1000 Touch thermal cycler.

BIO-RAD

5. If the version installed on your instrument does not match the current version listed on the website, download and install the update.

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.

Note: CFX Manager software version 2.1 can automatically check for updates if the computer is connected to the Internet. For more information, refer to the CFX96 Touch™ and CFX384 Touch™ Real-Time PCR Detection Systems Instruction Manual.

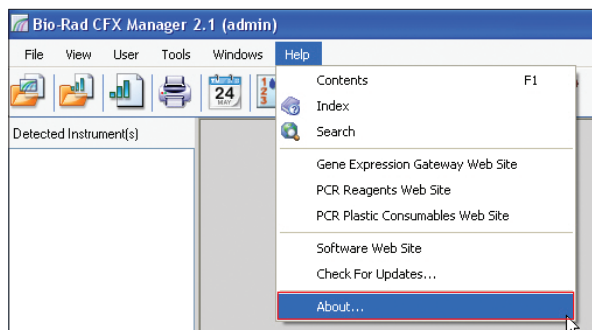


Fig. 4. Navigating in the main window of CFX Manager software.

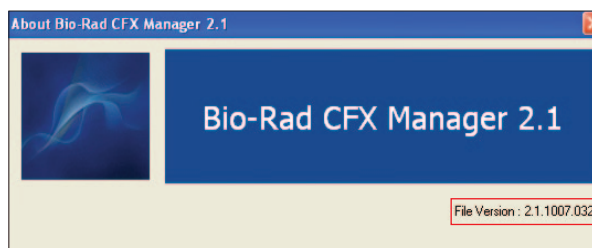


Fig. 5. Installed software version displayed in the About Bio-Rad CFX Manager 2.1 window.

Notice regarding Bio-Rad thermal cyclers and real-time systems:

Purchase of this instrument conveys a limited non-transferable immunity from suit for the purchaser's own internal research and development and for use in human in vitro diagnostics and all other applied fields under U.S. Patent Number 5,475,610 (Claims 1, 44, 158, 160–163, and 167 only), or corresponding claims in its non-U.S. counterpart, owned by Applied Biosystems. No right is conveyed expressly, by implication, or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applied Biosystems' U.S. Patent Number 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 Numbers 6,767,512 and 7,074,367.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 31 3689 6600
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Malaysia** 60 3 2117 5260 **Mexico** 52 555 488 7670
The Netherlands 0318 540666 **New Zealand** 64 9 415 2280 **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 3170 **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 **Thailand** 66 2 6518311 **United Kingdom** 020 8328 2000