



# CFX Opus 96, CFX Opus 384, and CFX Opus Deepwell Real-Time PCR Systems

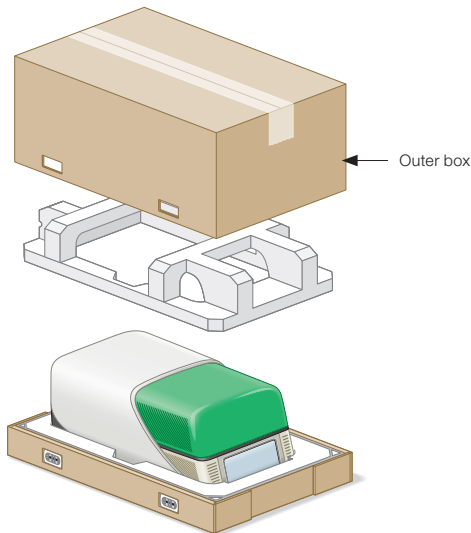
## Quick Start Guide

### Unpacking and Installing the CFX Opus System

The CFX Opus System is packed in a single box with the accessories included in the box. The accessories kit includes the power cord, USB cable, shipping screw plug, and an Ethernet cable. Set the box on the benchtop and follow the instructions per the labels affixed to the shipping box. For more detailed instructions, please refer to the instrument guide.

#### Unpacking the CFX Opus System

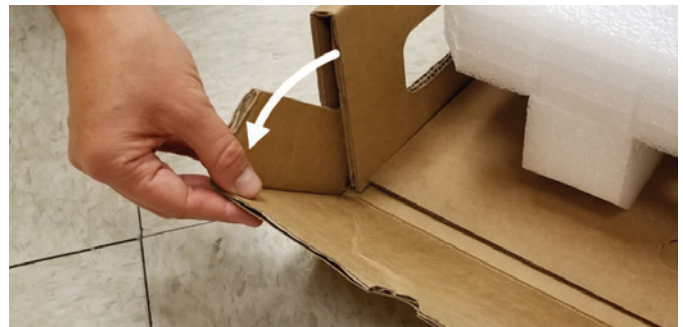
1. Remove the straps from the shipping box.  
DO NOT cut the top of the box open. Go to step 2 to continue unpacking the instrument.



2. Pinch and pull out the clips. After pulling out the four clips, lift up the outer box to expose the instrument.



3. Pull down for tray access.

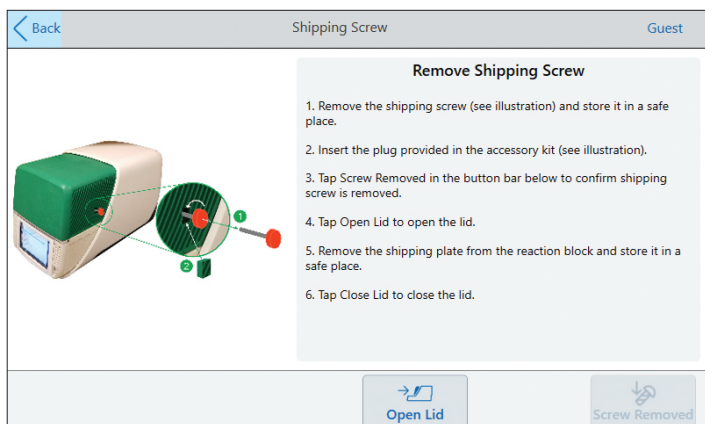


4. Slide tray to pull out the instrument.



### Removal of the Shipping Screw

1. Use the supplied power cord to plug the CFX Opus System into an appropriate electrical outlet.
2. Turn on the CFX Opus System using the power switch located on the right-hand side on the back of the instrument.
3. Upon power-up, follow the instructions displayed on the touch screen to remove the shipping screw, which is located on the right-hand side of the instrument.



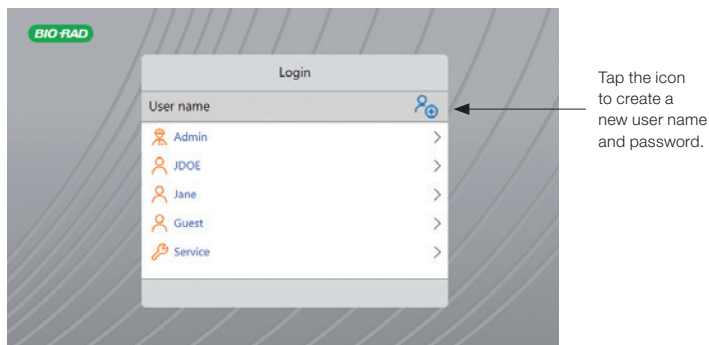
### Overview of CFX Opus Software

For additional details about CFX Opus Software, please refer to the instrument guide (10000119983).

### Login Screen

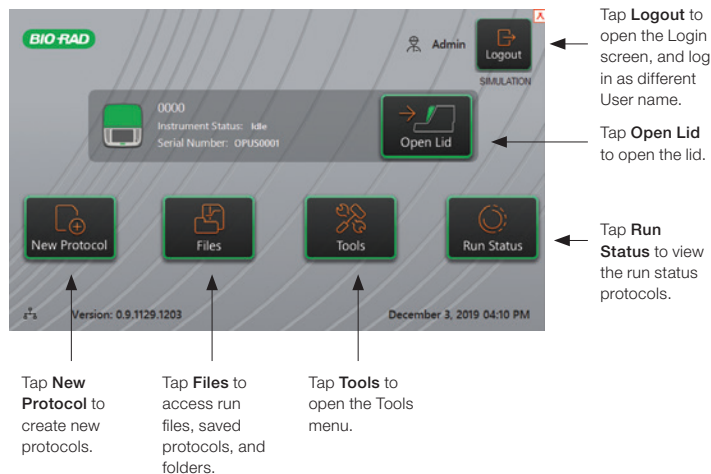
Create new user names by tapping the icon, or log in as Admin or Guest.

The Service login user name is reserved for Bio-Rad Service personnel.



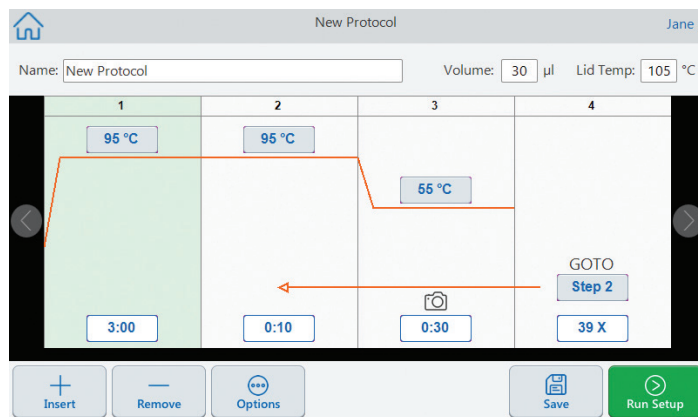
### Home Screen

Tap the buttons in the Home screen to perform the functions and to view connectivity and instrument status.



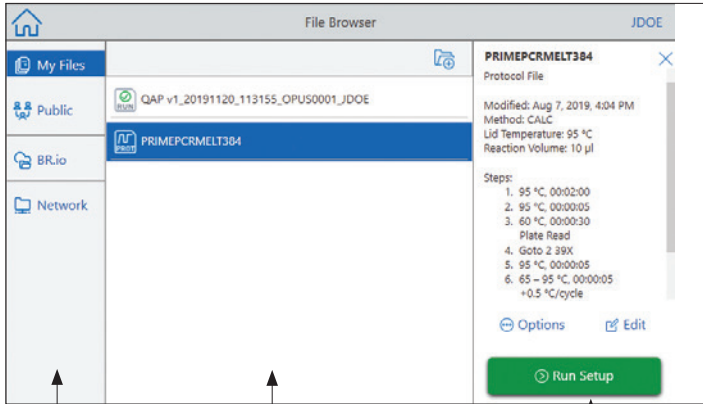
### New Protocol

Tap the buttons in the bottom bar in New Protocol to insert, remove, or modify steps in the protocol window. Tap **Save** to save any changes to the protocol. The camera icon indicates a "Read." To adjust, tap **Options**. Tap **Run Setup** to continue the process of starting a run.



**Files**

Tap **Files** in the Home screen to open File Browser. Tapping on a directory on the left-hand side, for example, MyFiles, will display the list of files (protocols, completed runs) in the center of the File Browser screen. Tapping on a specific file type will display the metadata information on the right-hand side of the File Browser screen. From the metadata pane, you can copy or delete a file from Options, or edit a protocol by tapping **Edit**.



Directory list

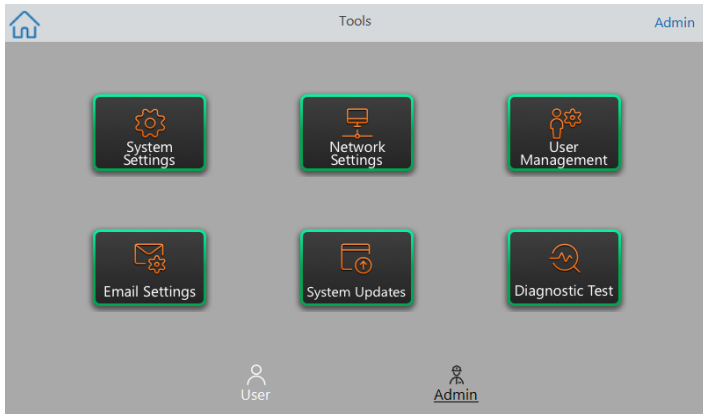
File list

Metadata pane

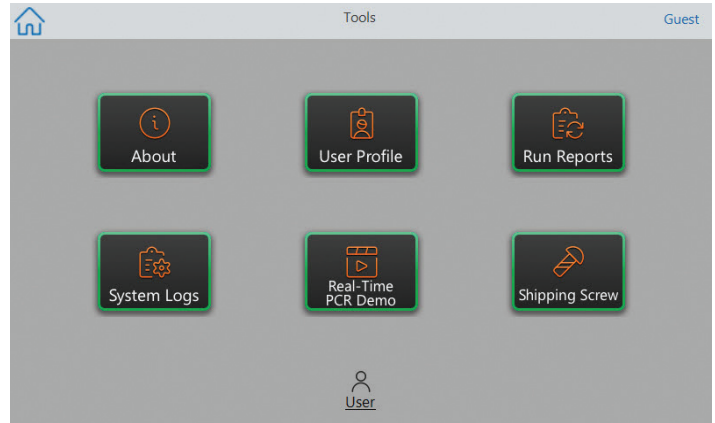
**Tools**

Tap **Tools** in the Home screen to open the Tools menu. There are two possible displays depending on whether you are logged in as Admin or as a User.

As an Admin, you are able to access the Admin and User screens as shown.



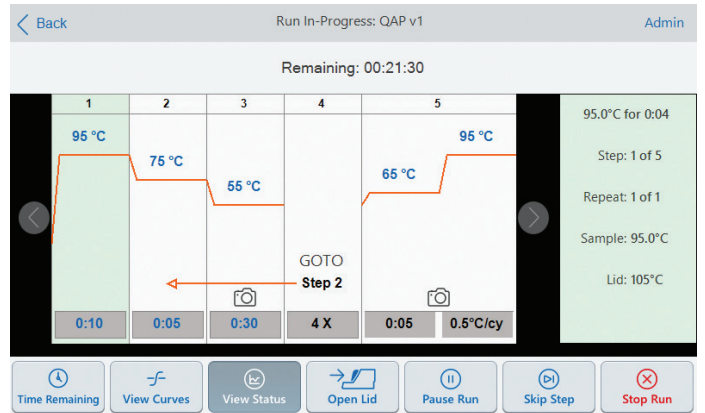
Logged in as Admin, you have access to the Admin tools.



Logged in as Guest or User, you have access to the User tools.

**Run Status**

Tap **Run Status** in the Home screen to open Run Status. You can access and view your run in progress by tapping the buttons in the bottom bar.



For more details, refer to the instrument guide (10000119983).

Visit [bio-rad.com/CFXOpus](https://www.bio-rad.com/CFXOpus) for more information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc.  
All trademarks used herein are the property of their respective owner.



**Bio-Rad  
Laboratories, Inc.**

---

Life Science  
Group

**Website** [bio-rad.com](https://www.bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23  
**France** 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050  
**Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670  
**The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23  
**Russian Federation** 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23  
**Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

