
Upgrading ChemiDoc or ChemiDoc MP (2017, touch screen model) Imaging System Running Image Lab Touch version 2.0, 2.1, 2.2, or 2.3 to 2.4

Before You Begin

1. Obtain a USB drive and insert it into a USB port on the computer.
Tip: Bio-Rad recommends the FAT32 format.
2. Copy the file ImageLabTouchSetup.deb to the root directory on the USB drive.
3. Safely remove the USB drive from the computer.

Installing Image Lab Touch v2.4

1. Power on the ChemiDoc or ChemiDoc MP instrument.
2. When Image Lab Touch software starts, log in as the Admin user if one was created.
3. Insert the USB drive into a USB port on the instrument.
4. If the software version on the USB drive is higher than the software on the instrument, a message appears prompting you to upgrade the software.

Touch OK to proceed with the upgrade.

Note: If the message does not appear, tap Update Installation on the System Settings menu on the right side of the toolbar. A dialog box appears displaying the current version of the software on the instrument and the update version on the USB drive. If the version on the USB drive is newer, proceed with the upgrade.

5. To complete the installation, follow the instructions that appear on the touch screen.
When the installation completes, you can remove the USB drive from the instrument.

Contacting Technical Support

Bio-Rad Technical Support in the United States is open Monday–Friday, 5:00 AM to 5:00 PM, Pacific Time.

Phone: 1-800-424-6723, option 2

E-mail: Support@bio-rad.com (U.S./Canada Only)

For technical assistance outside the U.S. and Canada, contact your local technical support office or click the Contact us link at www.bio-rad.com.

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions.

All trademarks used herein are the property of their respective owner.

Copyright © 2018, Bio-Rad Laboratories, Inc. All Rights Reserved.