
Upgrading ChemiDoc and ChemiDoc MP Imaging Systems Running Image Lab Touch Software 2.3 or 2.4 to Image Lab Touch Software 3.0.1

New Features in Image Lab Touch Software 3.0.1

Image Lab Touch Software 3.0.1 includes the following new features:

- Support for Simplified Chinese language
- Ability to print or export an image using the current display zoom area
- Simplified user interface for entering the network path to export files
- Support for future feature enhancements
- Increased onboard storage of 15–20 GB for gallery images

Upgrading Image Lab Touch Software to Version 3.0.1

IMPORTANT: After the Imaging System is upgraded to Image Lab Touch Software version 3.0.1, it cannot be downgraded to an earlier version.

Tip: Although all user data and images will remain on the imaging system after the upgrade, Bio-Rad recommends that you back up all gallery images by exporting them to a separate USB drive. Bio-Rad also recommends that you remove unwanted images from the imaging system before upgrading to shorten the upgrade time.

1. Power on the ChemiDoc or ChemiDoc MP Imaging System.
2. When Image Lab Touch Software starts, log in as the Admin user if one was created.
3. Insert a USB drive with sufficient free space to export all gallery images into a USB port on the Imaging System and export all gallery images.

IMPORTANT: Do not export the images to the Image Lab Touch Software v3.0.1 USB drive.

4. After the export completes, remove the USB drive.
5. Insert the Image Lab Touch Software v3.0.1 USB drive into the USB port on the imaging system's front panel.

6. Tap the Settings dropdown menu on the main toolbar and select Update Installation.
7. Follow the onscreen prompts to continue the upgrade.

Note: Depending on the number of images stored on the imaging system, the upgrade process can take up to 3 hours to complete. Do not power down or interrupt the imaging system during the upgrade.

After the upgrade is complete, the imaging system will automatically reboot. After the imaging system reboots, you can safely remove the USB drive.

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.

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

Copyright © 2020, Bio-Rad Laboratories, Inc. All Rights Reserved.