



Release Notes for ChemiDoc™ Touch Imager with Image Lab™ Touch Software

Version 1.2

February 2016

Contents

Contents.....	1
Introduction.....	1
New Features.....	2
Upgrading to Image Lab Touch 1.2 Software on the Imager.....	2
Documentation.....	3
Contacting Technical Support.....	4

Introduction

The ChemiDoc Touch imaging system is a stand-alone imaging instrument with Image Lab Touch software installed. The ChemiDoc Touch system supports acquisition of images from a wide range of laboratory gels and blots. It uses a sensitive camera with a charge-coupled device (CCD) and a large maximum aperture lens that provides high chemiluminescence sensitivity. The imager has a built-in UV transilluminator and white LEDs for epi illumination, and it works with UV, blue, and white trays that are used with gels and blots stained with a wide range of dyes and fluorophores.

The installed Image Lab Touch software enables you to acquire images in a three-step workflow: define the image size, select a gel or western blot application, and set the exposure. In addition, with Image Lab Touch, you can view images, fine-tune how images appear, print images, and export them to a computer running Image Lab software.

Images acquired with the ChemiDoc Touch imaging system can be exported and analyzed with Image Lab software.

New Features

Version 1.2 of the Image Lab Touch software includes these new features.

Camera zoom improvement

The ChemiDoc Touch camera zoom no longer returns to the fully zoomed-out position after you acquire an image, allowing you to acquire a new image with a different exposure time but the same image size. The ChemiDoc Touch camera zoom returns to the fully zoomed-out position only when you open the imager main door.

Ability to merge images

You can now merge two images of the same sample into a single image. A common example is combining a colorimetric image of prestained standards with a chemiluminescent image of the same blot. You can use the merged image to estimate the molecular weight or size of proteins on the chemiluminescent blot.

Upgrading to Image Lab Touch 1.2 Software on the Imager

Use this procedure to upgrade a previous version of Image Lab Touch installed on the ChemiDoc Touch imager to version 1.2.

1. Obtain a USB flash drive with the following characteristics:
 - Supports USB 2.0 or above
 - Is formatted with a recognized Windows format
 - Has no encryption software or other software add-ons on the drive
2. At the root of the flash drive directory structure, create a folder named Updater.
3. Log in to the Bio-Rad ChemiDoc Touch product web page and download the version 1.2 software zip file.
4. Extract the contents and copy the ImageLabTouchSetup.exe to the Updater folder on the flash drive.
5. Make sure the imager is turned on and Image Lab Touch is running.
6. Insert the USB flash drive into the USB port on the imager front panel. The system prompts you to confirm whether to proceed or not.
7. Press OK to proceed.

Caution: Do not interrupt the process. When the upgrade process is finished, the imager reboots automatically.

Documentation

The ChemiDoc Touch imaging system with Image Lab Touch software 1.2 release includes the following documentation, which has been updated:

- The ChemiDoc Touch Imaging System with Image Lab Touch Software User Guide, which is available on the Bio-Rad website ChemiDoc Touch product page. You must log in to access the documentation.
- Help, which is available on the imager touch screen.

The user guide for the ChemiDoc Touch imager with Image Lab Touch software 1.2 includes the following changes:

- Added explanation of the merge images feature plus instructions on how to merge two images
- Added instructions about how to export to an external hard drive and a shared folder on a stand-alone computer in the Exporting Images chapter
- Added Appendix C, Exporting to a Shared Folder, which explains how to set up a shared folder on a stand-alone computer
- Removed information about Image Lab software

The ChemiDoc Touch imager user guide contains information about the instrument and the installed Image Lab Touch software. Information about how to use Image Lab software to analyze images acquired with the ChemiDoc Touch system will be available in a new Image Lab software user guide to be released with Image Lab software 6.0. To read about Image Lab software, access the 1.0 version of the user guide for the ChemiDoc Touch imager, which is available on the Image Lab software version 5.2 Help menu.

Adobe Reader is required to open the user guide. If you do not already have Adobe Reader, go to acrobat.adobe.com/us/en/products/pdf-reader.html to download it for free.

How to access the user guides in Image Lab software

- In the Help menu, select User Guides, then select an imager from the list.

Contacting Technical Support

The Technical Support department in the United States is open Monday through Friday, 5:00 AM to 5:00 PM, Pacific time. Go to consult.bio-rad.com for worldwide technical support.

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

Web: consult.bio-rad.com

Email: Support@bio-rad.com (United States/Canada only)

For technical assistance outside the United States and Canada, contact your local technical support office.

Copyright © 2016, Bio-Rad Laboratories, Inc. All Rights Reserved. Adobe and Adobe Reader are trademarks of Adobe Systems Incorporated. Windows is a trademark of Microsoft Corporation.