



## Release Notes for Image Lab™ Touch Software 2.2, 2.1, and 2.0

August 2017

### Contents

|                                    |   |
|------------------------------------|---|
| Introduction.....                  | 1 |
| Improved Feature — 2.2 .....       | 1 |
| New Features — 2.1 .....           | 2 |
| New Features — 2.0 .....           | 3 |
| Documentation.....                 | 4 |
| Issues Resolved .....              | 5 |
| Contacting Technical Support ..... | 5 |

### Introduction

Image Lab Touch is a powerful, easy-to-use software package for image acquisition and basic image analysis of a wide range of laboratory gels and blots on supported Bio-Rad Laboratories imaging systems.

Image Lab Touch Software 2.2 is compatible with the ChemiDoc™ MP (2017 model), ChemiDoc (2017 model), and ChemiDoc Touch Imaging Systems. This version is pre-installed on ChemiDoc MP and ChemiDoc Imaging Systems.

ChemiDoc Touch Imaging Systems running Image Lab Touch 1.0 or 1.2 are compatible with 2.2. It is highly recommended that customers running 1.0 or 1.2 upgrade to 2.2.

### Improved Feature — 2.2

#### Stain Free Uniformity

Stain free uniformity has been improved in 2.2.

## New Features – 2.1

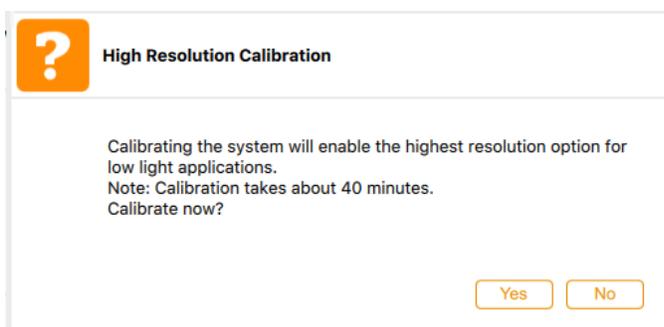
Image Lab Touch Software 2.1 adds the following new and improved features.

### New Setting for Chemiluminescent and Low-light Applications

A new highest resolution setting adds single-pixel resolution for chemiluminescent and low-light applications. Appendix A in the ChemiDoc and ChemiDoc MP Imaging Systems with Image Lab Touch Software User Guide includes detailed information about resolution and sensitivity.

### Highest Resolution Auto Exposure Calibration

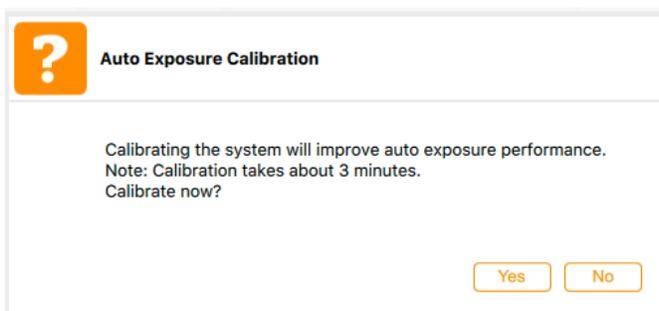
When the system starts up, it prompts the user to run a calibration. This calibration enables the new highest resolution setting for chemiluminescent and low-light applications.



The prompt appears each time the system starts up until the calibration is run.

### Auto Exposure Process Calibration

When the system starts up, it prompts the user to run a second, new calibration. This calibration enables Image Lab Touch 2.1 to detect too-faint signals more accurately.



The prompt appears each time the system starts up until the calibration is run.

## **Password Removal**

If no administrator has been assigned or if the administrator does not require Image Lab Touch users to log in with a password, users can remove their own password by entering a blank password in the Change Password dialog box.

## **New Features — 2.0**

Image Lab Touch Software 2.0 includes the following new and improved features.

### **Multiple User Accounts**

Each user can create their own user account where all their images are stored. Each user's last protocol used persists in the software. User accounts can be password protected by the user or by an administrator.

### **Advanced Administration Features**

The new administrator role is optional but strongly recommended. The administrator user can set parameters for other users by enforcing password protection, restricting exporting file options, and resetting user passwords. The administrator can also access the images stored in other user accounts. These features help ensure increased data integrity and security and help organizations comply with US FDA 21 CFR Part 11 regulations, which govern electronic records.

### **Enhanced User Interface**

Image Lab Touch 2.0 includes a new, more streamlined user interface, which features two main views, a Camera view and a Gallery view. Users start Image Lab in the Camera view, in which they can view in real time the sample to be imaged. Following the workflow in the left pane, the user can rapidly set parameters for imaging a gel or blot. The Gallery displays all images a user has saved. Images can be found by searching for their specific tags and dates.

### **Quick Help Screens**

To acquaint users with new features, several quick help screens appear throughout the software. These screens appear only once but can be easily recalled.

### **Multichannel Acquisition Support**

Users of the ChemiDoc MP Imaging System can now acquire up to three images of the same blot at the same time using different imaging parameters (applications). Users of the ChemiDoc and ChemiDoc Touch Imaging Systems can acquire up to two images (chemiluminescent and colorimetric). Multichannel images are automatically saved on the imaging system's computer hard drive as a single file that consists of the individual images and one composite image.

## Additional Applications

As in previous versions of Image Lab Touch, users select an application parameter based on the type of sample (gel or blot) and detection method used in the experiment (for example, chemiluminescence). Many more applications have been added to Image Lab Touch 2.0, including a wide range of Epi-light fluorescence applications (StarBright™ Blue 700, DyLight 800, Cy5, and the like) for use on the ChemiDoc MP only.

## Image Preview

Before acquiring a final image, users can preview the sample to evaluate the selected parameters. This preview image is a quick exposure taken at a lower resolution than the final image. In the preview image, users can also specify a region of interest on the blot (or gel) for automatic exposure.

## Intensity Values Check

Users can now rapidly determine the signal intensity and percentage of saturation at a single point of an image by tapping on the point. A box displays the parameters.

## Enhanced File Export Options

Images acquired are automatically saved to the imaging system's computer hard drive. Saved images can be exported in one of four formats: .scn (Image Lab), .tiff, .tiff (raw), or .jpg. Image files can be exported via USB drive, networked hard drive, or attached PC shared folder.

## Enhanced File Naming and Gallery Organization

Images acquired by Image Lab Touch are saved and stored in each user account according to the naming parameters (tags) set by the user.

## Documentation

Image Lab Touch includes an online help system. Tap the question mark icon in the main toolbar to display a relevant help topic. Tap Help Topics at the bottom of any help topic to view a list of all available topics.

The following user guides are available for download as PDFs on bio-rad.com. For ChemiDoc MP and ChemiDoc user guides, go to <http://www.bio-rad.com/en-us/category/chemidoc-imaging-systems>. The ChemiDoc Touch user guide can be found on bio-rad.com by searching for ChemiDoc Touch. User guides for Image Lab Software can be found at <http://bio-rad.com/en-us/product/image-lab-software>.

The following user guides can also be accessed on the Image Lab Software Help menu.

- ChemiDoc MP Imaging System with Image Lab Touch Software User Guide
- ChemiDoc Touch Imaging System with Image Lab Touch Software User Guide
- Image Lab Software User Guide Version 6.0

## Issues Resolved

The following issues are resolved in v. 2.2.

### **White Tray Calibration Could Impact LED Power Balance for Fluorescent Multiplex Applications**

Calibrating the white sample tray in Image Lab Touch 2.0 and 2.1 could impact the LED power balance. LED power balance is needed to provide the best data quality for fluorescence western blotting. This issue is resolved in 2.2. Impacted units will be corrected.

### **White Tray Calibration During Instrument Warmup Not Saved**

The system might not save white tray calibration performed during instrument warmup, making it necessary to repeat the calibration after warmup.

### **Renaming SAM Images Could Cause System Reset**

Occasionally, renaming a large number of images could cause the system to reset. This issue was seen most often when renaming SAM images.

### **Bright Pixels Appear in SAM Acquisition**

In SAM image acquisition bright pixel correction was not being applied, causing artificially bright pixel artifacts to appear in some images.

## Contacting Technical Support

The Bio-Rad Technical Support department is open Monday through Friday, 5:00 AM to 5:00 PM, Pacific time.

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

Email: [Support@Bio-Rad.com](mailto: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](http://www.bio-rad.com).

Cy is a trademark of GE Healthcare. DyLight is a trademark of Thermo Fisher Scientific Inc.

Copyright © 2017, Bio-Rad Laboratories, Inc. All Rights Reserved.