



## Release Notes for Image Lab™ Touch Software

### Version 2.0

February 2017

### Contents

Contents .....	1
Introduction.....	1
New Features .....	1
Documentation .....	3
Contacting Technical Support.....	4

### 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 Version 2.0 is compatible with the ChemiDoc™ MP (2017 model), ChemiDoc (2017 model), and ChemiDoc Touch Imaging Systems. This version is pre-installed on the ChemiDoc MP and ChemiDoc Imaging Systems.

ChemiDoc Touch Imaging Systems running Image Lab Touch v1.0 or v1.2 are compatible with v2.0. It is highly recommended that customers running these versions upgrade to v2.0.

### New Features

Version 2.0 of Image Lab Touch software includes the following new and improved features.

#### Multiple User Accounts

Each user can create their own user account where all their images are stored. Their last protocol used persists in the software. These 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 CFR 21 part 11, which governs electronic records.

### **Enhanced User Interface**

Image Lab Touch v. 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 gel or 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 v. 2.0, including a wide range of Epi-light fluorescence applications (StarBright 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 2.0 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, V. 6.0

## **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).

Adobe and Adobe Reader are trademarks of Adobe Systems Incorporated. DyLight is a trademark of Thermo Fisher Scientific Inc.

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