



Release Notes for Image Lab Touch Software 3.0.1 and 2.4

July 2020

Contents

Contents.....	1
Introduction.....	1
New Features — 3.0.1	2
Resolved Issues — 3.0.1	2
Improved Feature — 2.4	3
Documentation.....	3
Contacting Technical Support	3

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 3.0.1 is compatible with the 2017 model ChemiDoc and ChemiDoc MP Imaging Systems. This version is preinstalled on new ChemiDoc and ChemiDoc MP Imaging Systems. Customers running Image Lab Touch Software 2.3 or 2.4 on their existing ChemiDoc and ChemiDoc MP Imaging Systems can upgrade to 3.0.1.

Note: Image Lab Touch Software 3.0.1 is not compatible with GelDoc Go Imaging Systems, which ship with version 3.0 preinstalled.

New Features — 3.0.1

Support for Chinese (Simplified) Language

Image Lab Touch Software 3.0.1 now supports Chinese (Simplified) language. The software can be set to display the user interface in the Simplified Chinese language. It also provides online help in the Simplified Chinese language. Additionally, users can download the Image Lab Touch Software User Guide in Simplified Chinese from Bio-Rad's website.

Language-Specific Onscreen Keyboard

Image Lab Touch Software 3.0.1 can be set to display the onscreen touch keyboard in the Simplified Chinese language.

Improved Network Configuration User Interface

The user interface for setting the static IP address has been improved in Image Lab Touch Software 3.0.1.

Ability to Print the Current View

Users can now print or export the current display of an image (including the zoomed area) in Image Lab Touch Software 3.0.1.

Increased Internal Storage

After upgrading to Image Lab Touch Software 3.0.1, current ChemiDoc and ChemiDoc MP Imaging Systems will have approximately 20 GB of additional disk space.

Resolved Issues — 3.0.1

The following issues have been resolved in Image Lab Touch Software 3.0.1:

- On systems on which an Admin user has been defined, only those users with the Admin role can change the time or time zone.
- Image Lab Touch 3.0.1 supports all global time zones, resolving the issue that Daylight Saving Time settings in some locales were not supported.
- The ability to recover corrupted System Log files has been improved.

Improved Feature — 2.4

Image Lab Touch 2.4 allows ChemiDoc and ChemiDoc MP users to export to network servers and newer Windows 10 systems that have more restrictive security policies.

Note: Connecting to an SMB3 file server via a DFS namespace or alias will fail. You must use the fully qualified file server name. This is a known issue with Linux systems. Contact your IT department for assistance.

Documentation

Image Lab Touch Software 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 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 bio-rad.com/en-us/product/image-lab-software.

Users can also access the Image Lab Software User Guide version 6.1 from the Help menu in Image Lab Software.

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