



## Release Notes for Image Lab Software

### Version 6.0

March 2017

## Contents

Contents.....	1
System Requirements .....	2
New Features.....	2
Issues Resolved .....	3
Documentation.....	4
Bio-Rad Technical Support Department .....	4

## Introduction

Image Lab™ desktop software is a powerful application for image acquisition and analysis of a wide range of laboratory gels and blots on supported Bio-Rad Laboratories imaging systems. This version of the software also has advanced features for the normalization of western blots using either total protein normalization or housekeeping proteins. Image Lab 6.0 is free of charge for existing Bio-Rad ChemiDoc™, GelDoc™, and GS-900™ users.

Image Lab 6.0 is used for data analysis of images generated on the ChemiDoc MP, ChemiDoc, ChemiDoc XRS+, GelDoc XR+, and GelDoc EZ Imaging Systems along with the GS-900 Calibrated Densitometer. Images captured by the discontinued ChemiDoc MP (non-touch screen, 2011 model) and ChemiDoc Touch (2014 model) Imaging Systems can also be analyzed in Image Lab 6.0.

Image Lab 6.0 is also used for image acquisition on all non-touch screen Bio-Rad Imaging Systems (ChemiDoc XRS+, GelDoc XR+, GelDoc EZ, GS-900, and the discontinued ChemiDoc MP).

Image acquisition on all touch-screen Bio-Rad imaging systems is performed using Image Lab Touch 2.0, another member of the Image Lab software family.

10000076658 Rev A

## System Requirements

Install Image Lab 6.0 on a PC or Mac computer that meets the following requirements.

Component	Windows	Mac
Operating system	Windows 7 (32- and 64-bit) Windows 10 (64-bit)	OS X 10.11, 10.12
Processor	Intel Core i3 or equivalent	Intel Core i3 or equivalent
Hard disk space	60 GB	60 GB
Memory (RAM)	4 GB	4 GB
Screen resolution	1280 * 1024 or higher, 128 MB video RAM	120* 1024 or higher, 128 MB video RAM

Image Lab 6.0 does not support instrument control on ChemiDoc MP (non-touch screen, 2011 model), ChemiDoc XRS+, and GS-900 instruments on Mac operating systems. However, you can analyze images acquired on these imagers in Image Lab 6.0.

## New Features

Version 6.0 of Image Lab desktop software includes the following new and improved features.

### Increased Compatibility

Image Lab is now compatible with Windows 7 (32-bit and 64-bit), Windows 10 (64-bit), and Mac OS 10.11 and 10.12.

### Streamlined Lane and Band Editing

Lane and band editing has been streamlined by making lane editing more modeless. Users no longer have to switch between width, move, and bend modes. Users can now select and manipulate more than one lane at a time.

### Enhanced Normalization

Users can perform total protein normalization or housekeeping protein normalization more quickly and in fewer steps. It is now easier to link images for analysis. Any lane can now be defined as the normalization reference lane. Automated lane finding now detects the normalization channel and uses that lane frame for all the sample channels. Users can now copy lane changes from the normalization channel by pressing a button. A new lane statistics table, including total lane volume, can be exported to present analysis results outside of Image Lab.

## Lane Profile

As in the previous version, users can automatically identify lanes on a gel or blot image and then automatically detect bands. The Lane Profile tool is useful for fine-tuning the boundaries of bands in a lane and for subtracting background. Without leaving the Lane Profile window, users can now add and delete bands. They can also change and dynamically view changes in background removal using the new rolling disk slider.

## Issues Resolved

The following issues have been resolved.

<b>Issue Number</b>	<b>Description</b>
1511	Standard type selected by default for Nucleic Acid gels should be set to DNA-type standards
1698	Sort volume table by volume number rather than volume label for volumes with non-custom labels
2082	Image zoom lost after switching focus between images
2084	Unable to save the file after the file was moved to a different location using Finder
2119	Agarose gels should not be available for blot applications
2132	Export cropped image as PDF - overlays look terrible on the image in PDF
2459	No prompt to register with Bio-Rad in "upgrade software" scenario
2485	Selecting window as "active" does not restore window to view from "minimized"
2565	Image lab will crash if user cancels signing process for images acquired from signed protocol
2632	The edit menu, Select All" is enabled even when there aren't any items to select (Volume and Annotation Tools View)"
2633	The edit menu option, Paste" is always enabled even when there isn't a lane frame on the clipboard (Lane and Bands Tools)"
2694	Lane profile view is not restored as expected by user
2699	Unexpected "document signing cancelled" mean in an Image Lab Security Edition log?
2733	Image Lab crashes on Windows 10 Tech Preview
2736	Band can wrap and be displayed as negative number
2738	Unable to open .1sc files including unicode characters
2739	Can't rename permission groups or enable write once folders in non-English system
2740	Analysis table appears empty for image with Lanes/Bands detected when image is opened
2742	Adding a volume to image garbles log export output due to quotations messing up CSV fields.
2743	Customer attempts to "merge" images after splitting multichannel image and cannot do it (reported by Tech Support)
2747	Image lab using (um) instead of ( $\mu\text{m}$ ) as a symbol for micrometer units
2751	When creating a report for multichannel images, the band strip image from channel 1 is displayed for subsequent channels
2752	Band peak values change when band peaks recalculated after file re-opened

## Documentation

The following guides are available as PDFs in the Image Lab Software Help menu.

Image Lab Software User Guide, V. 6.0 — concepts and procedures for users of all Bio-Rad non-touch-screen imaging systems.

Information about each imaging system and how to start working with it now appears in an instrument guide:

- GelDoc EZ Imaging System with Image Lab Software Instrument Guide
- GelDoc XR+ and ChemiDoc XRS+ Imaging Systems with Image Lab Software Instrument Guide
- GS-900 Calibrated Densitometer with Image Lab Software Instrument Guide
- ChemiDoc MP (no touchscreen, discontinued) Imaging System with Image Lab Software Instrument Guide

Also included in the Help menu is ChemiDoc and ChemiDoc MP (touchscreen) Imaging Systems with Image Lab Touch Software User Guide

ChemiDoc and ChemiDoc MP touch screen instruments are controlled by Image Lab Touch software. Images obtained by Image Lab Touch can be exported to Image Lab 6.0 for more advanced analysis.

Up to date information on both Image Lab 6.0 and Image Lab Touch 2.0 can be found at <http://www.bio-rad.com/category/imaging-systems-software>

Adobe Reader is required to view the PDFs. To download a free version of Adobe Reader, go to <http://getadobe.com/reader/>.

## Bio-Rad Technical Support Department

The Bio-Rad 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](http://consult.bio-rad.com) for worldwide technical support.

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

**Web:** [consult.bio-rad.com](http://consult.bio-rad.com)

**Email:** [Support@Bio-Rad.com](mailto:Support@Bio-Rad.com) (United States/Canada Only)

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

Copyright © 2017, Bio-Rad Laboratories, Inc. All Rights Reserved. Adobe and Adobe Reader are trademarks of Adobe Systems Incorporated. Mac OS is a trademark of Apple, Inc. Windows is a trademark of Microsoft Corporation.