

IMAGING

Stain-Free Software Upgrade Quick Start Guide

Stain-Free Imaging on the Gel Doc™ XR+ and ChemiDoc™ XRS+ Systems

This guide will walk you through the steps to install Image Lab™ software version 5.1 beta.

Note: If you already have Image Lab software installed, then you will have to run the installer twice. First to uninstall the previous version, and then a second time to install version 5.1 beta. Alternatively, you may use Control Panel/Uninstall a Program to remove the previous version first.

1 Installing Image Lab software version 5.1 beta

- After you have downloaded Image Lab software, proceed to installing the software by double clicking **Setup**
- Click **Next** to launch the installation wizard
- Select **"I accept the terms"** and click **Next**
- Click **Next** to install Standard Edition

Note: If you are in a regulated environment, you may wish to install the Security Edition. You will need to have purchased a Security Edition license from Bio-Rad Laboratories, Inc. to install the Security Edition. If you already have the Security Edition and have activated your license, reactivation of Image Lab is not necessary.

- Click **Next** to install to the default location or change to install to a different folder
- Click **Install** to begin installation. Click **Finish** to complete installation. Installation of Image Lab software version 5.1 beta is now complete
- If you are upgrading a Gel Doc XR+ system, it is now stain-free enabled and ready to image stain-free gels. For upgrading Image Lab software on a ChemiDoc XRS+ system, a flat fielding calibration step will be required. Take a look at the camera on your imager if you are unsure which Bio-Rad imager you have (see Figure 1)



Gel Doc XR+ System ChemiDoc XRS+ System

Fig. 1. Gel Doc XR+ and ChemiDoc XRS+ camera locations.

2 Performing flat fielding calibration (optional step for Gel Doc XR+)

- To perform flat fielding you will need an orange fluorescent reference plate (catalog #170-8008)
Note: To obtain an orange fluorescent reference plate, required for flat fielding calibration, contact Bio-Rad tech support at: lsg.techserv.us@bio-rad.com or 1-800-4-BIORAD, option 2.
- Select **Edit** in the menu bar of Image Lab software
- Then select **Instrument Setup**
- In the Flat Field section of the Instrument Setup window click **Reset** (see Figure 2)
- Click **Skip** to bypass the lens flat fielding calibration because this procedure was done during installation of the imager

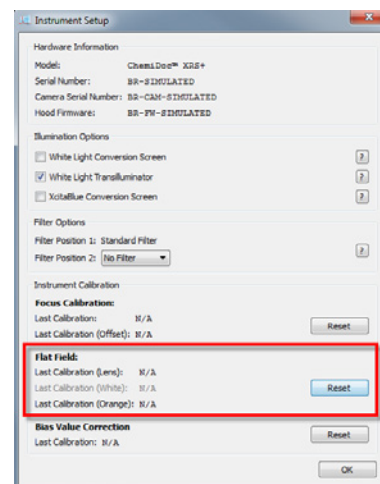


Fig. 2. Instrument Setup window.

- Go to the imaging hood and center the orange plate on the imaging stage, then close the door



- Click **Start Calibration**. On the upper right side of the hood, move the filter to position 1, then click **OK**. The flat fielding procedure will take a few minutes



- When the procedure is complete remove the orange plate. Click **OK** two times to exit out of the Instrument Setup window

Congratulations! Your Bio-Rad imager is now stain-free enabled. This additional calibration step generates flat field correction profiles needed to account for nonuniformity of the UV light source and is critical for accurate stain-free total protein normalization.



BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 11 5044 5699
Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00
France 01 47 95 69 65 Germany 089 31 884 0 Greece 30 210 9532 220 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300
Israel 03 963 6050 Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666
New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04
Singapore 65 6415 3188 South Africa 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 026 674 55 05
Taiwan 886 2 2578 7189 Thailand 800 88 22 88 United Kingdom 020 8328 2000