



IMAGING AND ANALYSIS

Getting Started with the GelDoc Go Imaging System

Quick Guide

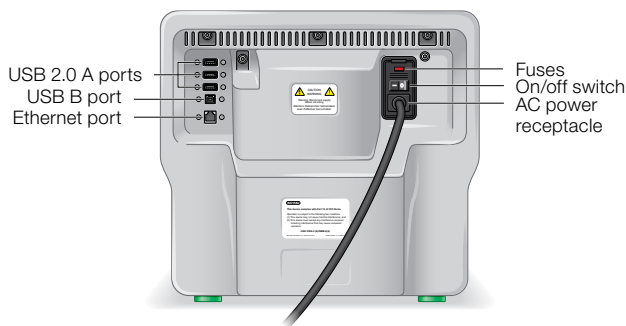
Set Up the Environment

The imaging system requires a space 38 x 37 x 46 cm (W x H x D) and a clearance of at least 8 cm from the back for instrument ventilation and for accessing the AC power cord. The imager should be placed on a sturdy and level laboratory bench or table away from excessive heat and moisture. The imager's operating temperature range is 10–28°C. The imager contains a universal power supply that supports a voltage range of 100–240 VAC.

Set Up the Imager

The imager ships with a power cable that works with AC 115/240 VAC outlets. The female end of the cable fits in the power receptacle located at the back of the imager.

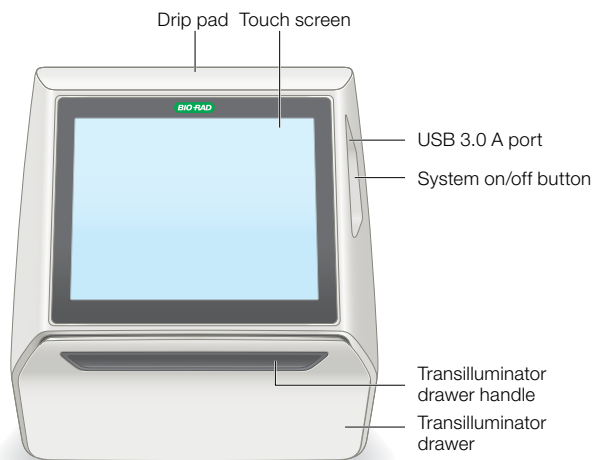
- 1 Locate the AC power receptacle on the back of your imager and connect the power cable to it. Plug the other end of the power cable into an AC outlet and turn the power switch to the **on** position.



- 2 For network access, connect an Ethernet LAN network cable to the Ethernet port on the back of the imager. Connect the other end of the cable to an Ethernet network port.

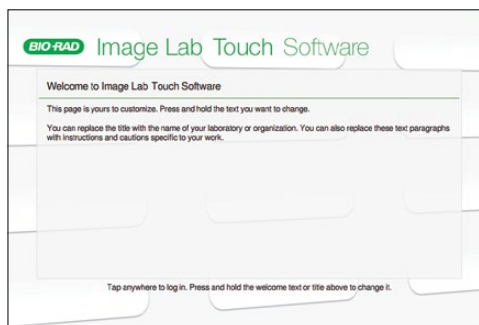
Note: When you save an image, you can specify the path to the network folder in which you want to save your image files. You can save multiple path names. Each saved path name appears in the dropdown list of destinations when you tap **Save to Network**.

- 3 On the front of the imager press the system on/off button to turn on the imager. Turning on the imager initializes the camera, computer, and Image Lab Touch Software.



First Time Login

When the imaging system is turned on, Image Lab Touch Software starts and the Welcome screen appears.



- 1 Tap anywhere on the touch screen.
- 2 Using the onscreen keyboard, tap in a user name and then tap **Create**.

The system prompts you to choose whether you want to be an administrator and displays information that helps you decide.

Note: An administrator must log in with a password. If no administrator has been specified, users can log in without a password.

- 3 If you do not want to be an administrator, tap **No**.
- 4 To become an administrator, tap **Yes** and create a password in the box that appears.

The Administrator Options dialog box appears, in which you can choose options for managing Image Lab Touch user accounts.

- 5 Choose the options you want and tap **Close**.

A Camera view of the imaging stage appears.

Check the Imager Setup

Check to ensure the following:

- The imager has been placed on a laboratory work bench surrounded by the required amount of space
- The imager is turned on
- Camera view is visible

Level the Imager

Is the imager level? In order to work properly, the imager must be situated on a level surface such as a laboratory work table. If the imaging bed in the imager is not level, samples — especially gel samples — might move during imaging.

To check whether the imager is level:

- 1 Locate the bubble level in the Imager Accessory Kit.
- 2 Pull the transilluminator drawer door to open.
- 3 Place a sample imaging tray on the imaging stage.
- 4 Place the bubble level at the approximate center of the sample tray.
- 5 If the entire bubble is visible inside the level's center ring, the imager is level and no adjustments are necessary.



- 6 If the entire bubble is not visible inside the level's center ring, adjust the level by rotating the leveling feet of the instrument to raise or lower it until the bubble appears at the center of the level.
- 7 Check whether the entire bubble is visible inside the level's center ring.
- 8 When the entire bubble is visible inside the center ring, remove the level from the sample tray.

Set Up the Printer

An optional Mitsubishi Thermal Printer (catalog #1708089) can be used with the imaging system.

To set up the printer:

- 1 Connect the printer's power cord to an electrical outlet.
- 2 Connect one end of the USB cable to the USB port on the back of the printer.
- 3 Connect the other end of the USB cable to one of the USB ports on the back of the imager.

Warning: Transilluminators are powerful sources of UV radiation, which can cause serious damage to unprotected eyes and skin. The accessory UV Shield (#12012164) provides some UV protection. However, the shield does not protect others standing in the area around the imager. In this configuration, the system is rated as Risk Group 3 (most hazardous) for actinic UV (200–400 nm) per IEC 62471:2009 "Photobiological safety of lamps and lamp systems." When viewing around the UV shield at a distance of 20 cm, the measured actinic UV emission is approximately 4 W/m², with a permissible exposure time of about 7 seconds.

Image Analysis Software

Images acquired on the GelDoc Go Imaging System can be analyzed for lanes/bands and volume quantification using Image Lab Software.

Licenses and upgrades for the Image Lab Software (available for Windows and Mac operating systems) are included in the purchase price of the instrument. Registered users can log in to their account or register as new users to download a free copy of Image Lab from the Bio-Rad website. Image Lab Software is also available for purchase on a USB drive. Visit [bio-rad.com/imagelab](https://www.bio-rad.com/imagelab) to learn more about the Image Lab family of software.

Find Out More

Visit [bio-rad.com/geldocgo](https://www.bio-rad.com/geldocgo) to learn more about the instrument and to download the full version of the user guide.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00
Germany 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600
Japan 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30
Poland 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580
Sweden 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300
United Kingdom 44 01923 47 1301

