

The Bio-Rad Imaging System Family

Unmatched Performance, Ease of Use, and Reliability for More Than 20 Years





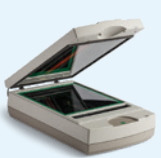
Bio-Rad provides a variety of imaging systems suitable for a wide range of applications. This guide can help you select the best system to detect and quantitate fluorescent, chemiluminescent, Stain-Free, and colorimetric blots and gels for your specific laboratory needs.

ChemiDoc Imaging Systems

Best-in-class performance with ease of use for fluorescence and chemiluminescence detection and all general gel documentation applications.

Gel Imaging Systems

Automated gel imaging and documentation for research laboratories and current good manufacturing practice environments.

	ChemiDoc MP	ChemiDoc	ChemiDoc XRS+	GelDoc Go	GS-900
					
	High-end imaging system for the best fluorescence and chemiluminescence detection	Standard system with chemiluminescence sensitivity matching X-ray film	Reliable, easy-to-use system with good chemiluminescence sensitivity	System for gel documentation in three simple taps for gels and colorimetric blots	High-resolution scanner for stained gels and blots
Blot Detection					
Chemiluminescence	Best	Best	Good	No	No
Fluorescent blots	Full RGB, far red, near infrared	Blue only	No	No	No
Multiplex fluorescence	Up to 3 dyes	Upgradeable	No	No	No
SYPRO Ruby Protein Blot Stain*	Yes	Yes	No	Yes	No
Colorimetric blots	Yes	Yes	Yes	Yes	Yes
Stain-Free blots**	Yes	Yes	Yes	Yes	No
Nucleic Acid Detection					
Ethidium bromide	Yes	Yes	Yes	Yes	No
SYBR®, GelGreen, and GelRed Stains	Yes	Yes	Yes	Yes	No
Fast Blast DNA Stain	Yes	Yes	Yes	Yes	Yes
Protein Gel Staining					
Stain-Free gels**	Yes	Yes	Yes	Yes	No
Coomassie Brilliant Blue, Coomassie Fluor Orange, and Silver Stains	Yes	Yes	Yes	Yes	Yes
SYPRO Ruby, Flamingo, and Oriole Gel Stains	Yes	Yes	Yes	Yes	No
Pro-Q Stain	Yes	Yes	Yes	No	No
Cy2, Cy3, Cy5 labels	Yes	Upgradeable	No	No	No
Specifications					
User interface	12 in. graphical touch screen	12 in. graphical touch screen	Keypad entry or computer control	9.7 in. multitouch display	Computer control
Maximum image area (W x H)	21 x 16.8 cm	21 x 16.8 cm	25 x 26 cm	21 x 14 cm	29 x 33 cm
Detector	Cooled CCD, 6 megapixels	Cooled CCD, 6 megapixels	Supercooled CCD, 1.4 megapixels	CMOS, 6.3 megapixels	Trilinear CCD
Dynamic range	>4 orders of magnitude	>4 orders of magnitude	>4 orders of magnitude	>3.5 orders of magnitude	0–3.4 OD transmissive absorbance
Smart Tray Technology	Yes	Yes	No	Yes	No
Auto-focus and auto-exposure	Yes	Yes	Yes	Yes	No
IQ/OQ kit	Yes	Yes	No	Yes	Yes

* Optimal with low fluorescence polyvinylidene difluoride membrane.

** Requires Bio-Rad Stain-Free Gels.

CCD, charge-coupled device; CMOS, complementary metal oxide semiconductor; IQ, installation qualification; OD, optical density; OQ, operational qualification; RGB, red, blue, and green.

Visit bio-rad.com/ImagingSystems to learn more about the Bio-Rad family of GelDoc Go and ChemiDoc Imaging Systems.



Bio-Rad Imagers Powered by Image Lab Software

Powerful Acquisition and Efficient Analysis

Image Lab Software allows you to acquire and analyze publishable gel and blot images with ease. The robust analysis package provides automated detection of lanes and bands and a simplified normalization process with complete report generation.

Features:

- Convenient preview mode
- Auto-exposure for optimal image quality
- Intuitive lane/band analysis
- Publication-quality export of TIFF or JPEG files
- Total protein normalization
- Accurate molecular weight analysis
- Optimal light source that can be selected
- Manual and automated calculation of exposure time

Imaging System	Image Lab Touch Software	Image Lab Software for Windows	Image Lab Software for Mac	Image Lab Software, Security Edition, for Windows*
ChemiDoc MP	Onboard acquisition	Desktop analysis	Desktop analysis	Desktop analysis
ChemiDoc	Onboard acquisition	Desktop analysis	Desktop analysis	Desktop analysis
ChemiDoc XRS+	No	Acquisition and analysis	Desktop analysis	Acquisition and analysis
GelDoc Go	Onboard acquisition	Desktop analysis	Desktop analysis	Desktop analysis
GS-900 Densitometer	No	Acquisition and analysis	Desktop analysis	Acquisition and analysis

* Researchers requiring a software package to help enable U.S. FDA 21 CFR Part 11 compliance can purchase Image Lab Software, Security Edition.

Ordering Information

Catalog # Description

ChemiDoc Imaging Systems

12003154	ChemiDoc MP Imaging System
12003153	ChemiDoc Imaging System
12003263	ChemiDoc to ChemiDoc MP Upgrade Kit
12005137	ChemiDoc and ChemiDoc MP Imaging System Installation/Operational Qualification Kit
1708265	ChemiDoc XRS+ Imaging System with Image Lab Software

GelDoc Go and GS-900 Systems

12009077	GelDoc Go Gel Imaging System with Image Lab Software
12012147	GelDoc Go Gel Imaging System IQ/OQ Kit
1707991	GS-900 Calibrated Densitometry System
1707994	GS-900 Installation/Operational Qualification Kit

Software

12012931	Image Lab Software on USB Drive, stand-alone version for PC or Mac
17006130	Image Lab Software, Security Edition, 1 license, U.S. FDA 21 CFR Part 11 module

Visit bio-rad.com/ImageLab to learn more about what Image Lab Software can do for you.

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



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399
Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23
France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300
Israel 0 3 9636050 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Luxembourg 00 800 00 24 67 23
Mexico 52 555 488 7670 The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23
Portugal 00 800 00 24 67 23 Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23
Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23 Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311
United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

