







The Bio-Rad Imaging System Family

Unmatched performance, ease of use, and reliability for over 20 years

At Bio-Rad, we are committed to providing a variety of cutting edge imaging systems that suit your needs for a wide range of applications. Detect and quantitate fluorescence, chemiluminescence, stain-free and colorimetric blots and gels with our systems, featured below.

	Gel Doc™ EZ	Gel Doc XR+	GS-900™	ChemiDoc™ XRS+	ChemiDoc	ChemiDoc MP
						
	Analyzes protein and DNA gels with the push of one button.	Analyzes larger protein and DNA gels.	High-resolution scanner for stained gels and blots.	Affordable system for chemiluminescent blotting and routine gel imaging.	Standard system for unmatched chemiluminescent performance and routine gel imaging.	The best-in-class all-in-one system for chemiluminescent and fluorescent blot imaging and routine gel imaging.
DNA gels (EtBr, SYBR®, GelRed, and more)	✓	✓	—	✓	✓	✓
Protein gels (Coommassie, stain-free, and more)	✓	✓	✓	✓	✓	✓
Fluorescent blots*	—	—	—	—	✓	✓
Multiplex fluorescent gels and blots*	—	—	—	—	—	✓
2-D protein gels	✓	✓	✓	✓	✓	✓
Colony counting	✓	✓	✓	✓	✓	✓
Chemiluminescent blots	—	—	—	✓	✓	✓
Colorimetric blots (Ponceau, AP, HRP)	—	—	✓	✓	✓	✓
Stain-free gels and blots	✓	✓	—	✓	✓	✓

* ChemiDoc MP Imager supports 400–800 nm range of all commonly used RGB and IR fluorophores (e.g., DyLight, Alexa Fluor, Cy, IRDye, and StarBright™ Blue 520 and Blue 700). ChemiDoc Imaging System is limited to blue-excitable fluorophores (e.g., StarBright Blue 520).

Visit bio-rad.com/imagingsystems to learn more about the Bio-Rad family of Gel Doc and ChemiDoc imaging systems.



Bio-Rad Imagers Powered by Image Lab™ Software

Powerful acquisition and efficient analysis

Image Lab Software is a frequently updated, user friendly, powerful software package for acquisition and analysis of gel and blot images.

Features:

- Free to download
- Convenient preview mode
- Auto exposure for optimal image quality
- Accurate molecular weight analysis
- Intuitive lane/band analysis
- Publication-quality export of TIFF or JPEG files
- Total protein normalization

Image Lab Touch and Image Lab Software run the Bio-Rad imaging system family.

Imaging System	Image Lab Touch Software	Image Lab Software* (PC and Mac compatible)	Compatible with U.S. FDA CFR 21 Part 11 Regulation
Gel Doc EZ Imager	—	✓ Acquisition and analysis	✓ Yes
Gel Doc XR+ Imager	—		
GS-900 Densitometer (PC only)	—		
ChemiDoc XRS+ Imager	—	✓ Analysis only	
ChemiDoc Imager	✓ Acquisition only		
ChemiDoc MP Imager			

* Image Lab Software includes analysis tools for normalization via total protein or your housekeeping protein.

Ordering Information

Catalog #	Description
1708270	Gel Doc EZ Imaging System
1708195	Molecular Imager® Gel Doc XR+ System
1707991	GS-900 Calibrated Densitometry System
1708265	ChemiDoc XRS+ Imaging System
12003153	ChemiDoc Imaging System
12003154	ChemiDoc MP Imaging System

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

Alexa Fluor, Dylight, and SYBR are trademarks of Thermo Fisher Scientific. Cy is a trademark of GE Healthcare. GelRed is a trademark of Biotium, Inc. IRDye is a trademark of LI-COR Biosciences. Mac is a trademark of Apple, Inc.



**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

