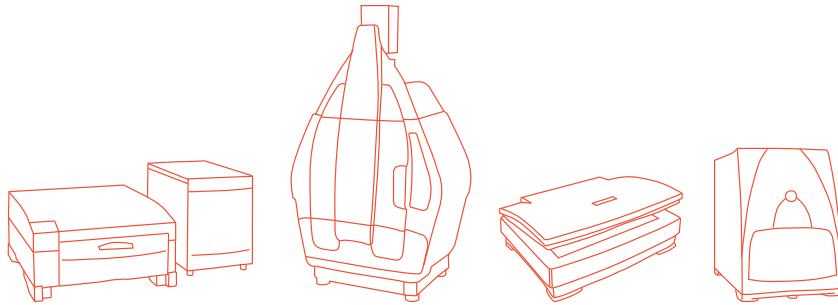




Bio-Rad Imaging Systems Family







Unlike other imagers, Bio-Rad imaging systems were developed by scientists for scientists. Bio-Rad understands your challenges and knows how to overcome them. That's why Bio-Rad imaging systems have consistently won the prestigious Life Science Industry Awards for most outstanding image analysis system. Life scientists choose the most reliable and best performing products to receive these awards in categories spanning all of the life science disciplines.

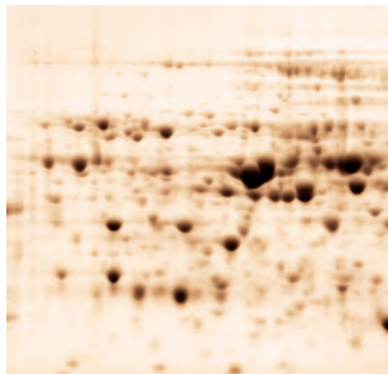


Bio-Rad Laboratories, Inc.

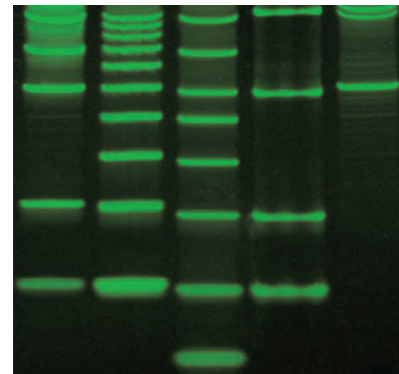
2006, 2007, and 2008 Award Winner
Best Image Analysis Systems

2010 Award Winner
Top Ten Innovations

COMPACT



EFFICIENT



DNA Gels



Protein Gels



3-Color Fluorescence



2-D Protein Gels



X-Ray Film



Colony Counting



2-D DIGE



Phosphor Imaging



Chemiluminescent Blots



Colorimetric Blot Detection



Stain-free imaging

Gel Doc™ EZ System



Imaging and analysis of protein and DNA gels in seconds with the push of just one button

About the system

- A simple and compact system that is customizable, easy to use, and that can fit easily on a laboratory bench
- Offers remarkable flexibility with its four tray options for different applications
- Provides expedited workflow and increased free time. A 2-hour Coomassie staining protocol can be condensed to 5 minutes using the Stain-Free tray on the Gel Doc EZ system
- Provides high-quality, publication-ready images without any manipulation or user-introduced errors
- Possesses the sophisticated and intuitive Image Lab™ software, which provides images and analyzed results that are highly reproducible, within a matter of seconds

Gel Doc™ XR+ System

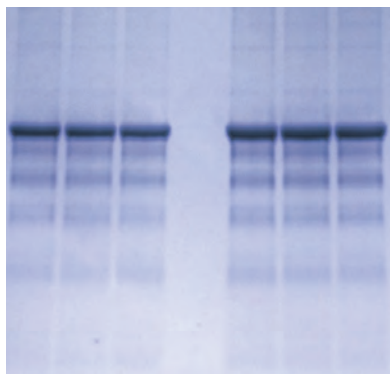


Affordable system for routine imaging of DNA and protein gels

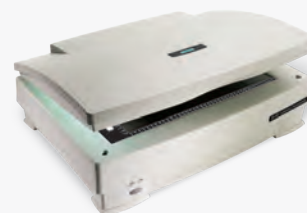
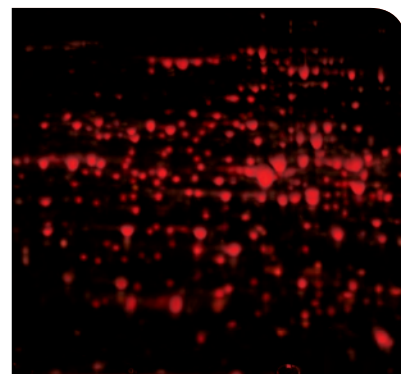
About the system

- Image Lab software provides automated and reproducible results from image acquisition to analysis
- Autofocus at any zoom level ensures high-quality images that are highly reproducible
- Trans-blue light plate allows on-plate band excision using safe DNA staining (optional)
- Can be upgraded to the ChemiDoc™ XRS+ system
- Installation qualification/operational qualification (IQ/OQ) protocols (optional)
- Quantity One® 1-D analysis software 21 CFR Part 11 module (optional)

AUTOMATED



POWERFUL



ChemiDoc™ XRS+ System



Affordable system for chemiluminescent western blotting and routine gel imaging

About the system

- System delivers both high sensitivity and excellent resolution
- Image Lab software provides automated and reproducible results from image acquisition to analysis
- Autofocus at any zoom level ensures high-quality images that are highly reproducible
- Trans-blue light plate allows on-plate band excision using safe DNA staining (optional)
- IQ/OQ (optional)
- Quantity One 1-D analysis software module is available for compliance with U.S. FDA 21 CFR Part 11 requirements (optional)

ChemiDoc™ MP System



Designed for chemiluminescence and multiplex fluorescence detection, this imager can also be used for routine gel imaging

About the system

- Our most sensitive charge-coupled device (CCD)-based imaging system
- Image Lab software provides automated and reproducible results from image acquisition to analysis
- Autofocus at any zoom level ensures high-quality images that are highly reproducible
- 3-color multiplex capability
- Automated, preprogrammable filter and light operation determined by the application selected
- Trans-blue light plate allows on-plate band excision using safe DNA staining (optional)
- Quantity One 1-D analysis software 21 CFR Part 11 module (optional)
- Total protein normalization with stain-free technology

GS-800™

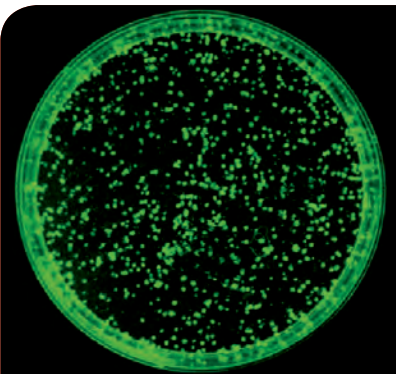
Calibrated Densitometer



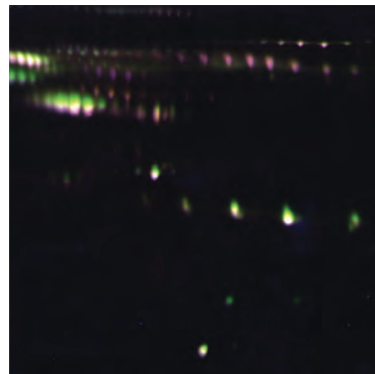
Affordable, high-resolution scanner for stained gels and blots

About the system

- Self-calibrating with internal standards for run-to-run consistency
- IQ/OQ (optional)
- Quantity One 1-D analysis software 21 CFR Part 11 module (optional)



REPRODUCIBLE



FAST



PharosFX™ Plus System



Laser-scanning system allows high-resolution 3-color imaging and phosphor imaging

About the system

- A high-resolution laser scanning system for fluorescent and radioisotopic applications
- 3-color multiplex capability
- Quantity One 1-D analysis software 21 CFR Part 11 module (optional)
- Fully upgradable from single laser to 3-laser capability

Personal Molecular Imager™ (PMI™) System



Basic phosphor imaging system for radioisotopic applications

About the system

- A high-resolution laser scanning system for radioisotopic applications
- Used with storage phosphor screens for ^{32}P , ^{33}P , ^{35}S , ^{14}C , and ^3H
- Quantity One 1-D analysis software 21 CFR Part 11 module (optional)

Imaging System Packages

For most researchers, no single imager addresses every experimental need. Therefore, Bio-Rad has created unified image acquisition and analysis software to allow one computer to operate a combination of instruments.

These are our most often requested imaging system combinations.

Gel, Blot, and Isotopic Imaging Package



▪ ChemiDoc XRS+ system

▪ Personal Molecular Imager (PMI) system

Multiplex Blot, Gel, and Isotopic Imaging Package



▪ ChemiDoc MP system

▪ Personal Molecular Imager (PMI) system

2-D Imaging Package



▪ PharosFX Plus system with blue and red lasers

▪ PDQuest™ Advanced 2-D analysis software

▪ EXQuest™ spot cutter

Life Science Industry Awards is a service mark of The Scientist.
Coomassie is a trademark of BASF Aktiengesellschaft.



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 31 3689 6600
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 30 210 777 4396 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 Malaysia 60 3 2117 5260 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 3170 South Africa 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700
Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 Thailand 66 2 6518311 United Kingdom 020 8328 2000