Imaging



Gel Doc™ XR+ Imager

Effortless and Accurate Gel Imaging and Analysis



Fast, Automated, Reproducible Imaging for Perfect Images Every Time

For more than two decades, the Molecular Imager[®] systems from Bio-Rad have been widely recognized and trusted high-quality imaging instruments. Whether you perform routine imaging of PCR products or protein purifications, the Gel Doc XR+ Imaging System will meet your needs. Designed for ease of use and an ability to support a wide array of applications, it is a perfect fit for individual laboratories, multiuser core facilities, or research on infectious diseases such as avian flu and H1N1 flu. Quality standards of engineering and manufacturing make the system adaptable to academic or biopharmaceutical laboratories. Thousands of researchers over the years have published Gel Doc XR+ System images in dozens of journals and established a record of reliability and excellence for the system.

The Gel Doc XR+ System enables quick and easy visualization, documentation, and analysis of nucleic acid and protein gels, blots, and macroarrays with a few clicks of the mouse. It is a flexible and easy to use system that supports fluorescence and colorimetric detection methods. The large imaging area accommodates a broad range of gel sizes, from large handcast polyacrylamide gels to small ReadyAgarose[™] Gels. The Gel Doc XR+ System is the ideal accompaniment to your PCR and electrophoresis systems, enabling image analysis and documentation of restriction digests, amplified nucleic acids, and genetic fingerprinting.

What the Gel Doc XR+ System Can Do for You:

Automated workflow from image to results

- Makes your gel imaging and analysis quick and effortless
- Eliminates the need for training
- Allows any user to repeat the same workflow in exactly the same way

BIO RAD

1

Gel Doc" XR+

Automated image capture

- Produces beautiful pictures of your gels
- Delivers quantitative analysis of your protein and DNA samples

Automated image optimization

 Delivers image data that is always optimized and reproducible without imaging artifacts

Automated image analysis

- Produces reports with data organized the way you want
- Minimizes time from lab work to presentations





SYBR[®] Safe Stain



Coomassie Brilliant Blue R-250 Stain



Oriole[™] Stain



Qdot blot

Gel Doc XR+ System Applications

Nucleic Acid Electrophoresis	Protein Gel Electrophoresis	Blotting
	Stain-free technology	
Ethidium bromide (EtBr)	Coomassie Blue	Colorimetric
SYBR [®] Green	Copper stain	Qdot 525
SYBR [®] Safe	Zinc stain	Qdot 565
SYBR [®] Gold	Flamingo [™] Fluorescent Gel Stain	Qdot 625
GelGreen	Oriole [™] Stain	Cy2
GelRed	Silver stain	Alexa Fluor 488
Fast Blast™ DNA Stain	Coomassie Fluor Orange	DyLight 488
	SYPRO Ruby	Fluorescein
	Krypton	



INS EP

Gel Doc XR+ System Is Powered with Image Lab[™] Software

Fast Results from a Completely Automated Workflow

The Gel Doc XR+ System is controlled by Image Lab Software to automatically and reproducibly generate gel images. Image Lab Software is fast, taking you from gel to printed results in seconds. Your results are visible with a single click of the mouse.

Image Lab Software eliminates the guesswork in imaging. You won't have to perform tedious repetitive steps to find the right focus setting. You won't have to guess which exposure times will best visualize the bands of interest. Image calculations and corrections are done automatically for your application. Whether you are working with protein gels, nucleic acid gels, or your own customized imaging application, the Gel Doc XR+ System will select the proper settings for optimum detection conditions of the stain or label.

You also won't need to set aside time for training new users; the system and the software are easy to use and work together from setup to finish. Imaging and image analysis can become the simplest work in the laboratory.

Simple and Reproducible Image Capture

Place your gel on the Gel Doc Imager's sample tray and run your protocol — your work is done in just one step. Design and save protocols for the imaging steps in your experiments and Image Lab Software will run the protocols exactly the same way every time. The protocol feature negates variability in results due to different people operating the imaging system. When a new project is related to an existing one, you can reuse the previously created protocol by revising its parameters and renaming it. Image Lab Software lets you devote your valuable time to research and discovery instead of to wondering if you used your imaging system properly.



Protocol Setup	1. Gel Imaging			
Acquisition Settings	Application			
	Nucleic Acid Gels Ethidium Bromide			
Analyze Image	Protein Gels			
2. Lane And Band Detection	Blots > SYBR Safe			
3. Analyze Molecular Weight	Custom SYBR Gold			
Report Settings	GelGreen			
4. Specify Content	GeRed			
	Select gel type: Bio-Rad Mini ReadyAgarose Gel ¥			
Protocol Summary	O Enter image area: 13.4 x 10.0 cm (WxiL)			
	Image Exposure			
	The software will automatically optimize the exposure time: Optimal Exposure Y			
	O Manually set exposure time: 0.500 sec			
Position Gel	Display Options			
Run Protocol	Highlight saturated pixels Image Color: Gray			

14

Automated workflow for any application.



The image stays focused at any zoom point.



Lane profiles depict band intensity and represent quantities of sample components separated in a gel.

Results and Reports

In addition to printing a picture of your gel for your records, Image Lab Software creates and prints reports of your experimental data. Any part of the report can be copied into popular document processing applications such as Adobe Acrobat and Microsoft Word, PowerPoint, or Excel files. To include a 3-D view of your gel, copy it using the Snapshot tool, and paste it into your presentation slide. High-quality, good-looking reports are easy to produce with the combined power of the Gel Doc XR+ System and Image Lab Software.

When an analysis parameter is changed, the results tables are updated instantly to reflect the new data.

Tutorials

With Image Lab Software, you don't need previous imaging experience to produce optimal gel images. Detailed tutorials are accessible via the toolbar and start-up page to acquaint you with all of the Image Lab Software capabilities.



Tutorials for novice users.

Sophisticated Software

Automated workflows — the entire workflow (image capture, results, report) is recorded in a protocol file. Protocols can be edited, resaved, reused, and shared among multiple users. Allows 100% repeatability of the workflow and ensures optimized image data and analysis specific to the selected application.

Auto focus — Image Lab Software's proprietary algorithms calibrate the system at setup for an automatic focused image at any zoom level. Eliminates user error and the need for manual camera adjustments to obtain an image, leading to higher image quality.

Auto camera aperture control — you do not have to focus or adjust aperture setting. Only adjust the zoom to position sample. Allows you to quickly image across different applications with different aperture setting requirements.

Flat fielding — flat fielding calibrations are performed for each application automatically. Delivers image data that is always optimized and reproducible without imaging artifacts for superior image uniformity and quantification.

Increased image resolution — decreases pixelation when images are cropped or zoomed. Yields smooth, clean images at any zoom level.



Images can be zoomed in without losing resolution.

Greater export functionality — no need to export images to another image editing program such as Photoshop Software to change the dpi before importing for publication. You can now define your desired resolution within Image Lab Software.

Molecular Biology Applications

The Gel Doc XR+ System can be used in multiple workflows. Gel electrophoresis remains an important tool for quantitative analysis of complex protein and nucleic acid samples. Applications include nucleic acid isolation or amplification for molecular cloning techniques, nucleic acid quality evalution prior to quantitative PCR, gene silencing, gene modulation, and gene expression. DNA gel electrophoresis is also used for fine separations of nucleic acid fragments or entire plasmids for analysis, cloning, transfection, environmental testing, and quality assurance testing. Gel Doc XR+ Systems are suitable for laboratories engaged in RNAi analysis, molecular diagnostics, epigenetics, pharmacogenomics, and forensic testing of genomic DNA.



E. coli colony screening by PCR.



ReadyAgarose 96 Plus Gel stained with EtBr.



PFGE gel stained with EtBr.

Safe DNA Imaging

Acquire images of DNA electrophoresis samples without exposing samples or users to harmful UV radiation by adding the XcitaBlue[™] Conversion Screen to your Gel Doc XR+ System. The screen enables detection of DNA with SYBR[®] Safe Stain and other blue light–excitable fluorescent stains such as GelGreen and SYBR[®] Green. Images and analyzed data produced by Gel Doc XR+ Systems are often incorporated into sample preparation and sample quantitation workflows.

Contraction line

Ethidium Bromide vs. SYBR® Safe DNA Stain Detection





Alternative to UV illumination to better preserve DNA samples. Top, serial dilutions of EtBr-stained precision molecular mass ruler (Bio-Rad Laboratories, Inc.) on agarose gel imaged with UV light; bottom, serially diluted precision molecular mass ruler stained with SYBR® Safe Stain on agarose gel imaged with XcitaBlue Conversion Screen. There is no loss in sensitivity when a combination of the nucleic acid SYBR® Safe Fluorescent Stain and less harmful blue excitation is used instead of UVexcitable EtBr. The image of the gel stained with SYBR® Safe Stain was taken using the XcitaBlue Conversion Screen and a filter for SYBR® Safe Stain and GFP. Less harmful detection methods preserve samples better and more efficiently for downstream uses such as cloning.

Gel Doc[™]XR+ Systems



Protein gels stained with silver stain (top) and Coomassie Blue Stain (bottom). Documentation of protein gels for lab notebooks or sample analysis, including protein purity and concentration assessment for differential protein expression studies or to monitor gene modulation products, is supported by the array of analysis tools in the Gel Doc XR+ System. The gel stained with silver stain shows salmon muscle, soybean, and rat brain extracts, protein mixtures, and *E. coli* extracts compared with Precision Plus Protein[™] Unstained Standards.

Protein gels fluorescence-stained with Oriole Stain (top) and Flamingo Fluorescent Gel Stain (bottom). Achieve higher levels of sensitivity with fluorescent stains to differentiate proteins with low expression. Resolve closely spaced spots or bands for protein profiling, quantitation, and characterization.

Capturing Gene Typing Results

Gel Doc XR+ Systems are the imagers of choice for pulsed field gel electrophoresis (PFGE). Gel Doc XR+ System images generated from CHEF Mapper[®] XA Systems for PFGE have been used often by PulseNet International (a global network of health and food regulatory agency laboratories coordinated by the Centers for Disease Control and Prevention) for strain typing and molecular epidemiology studies. Image data and analysis of PFGE gels have contributed to new findings in cancer research, food safety, public health, quality control, and genome mapping.



Images produced by the Gel Doc XR+ System have contributed to new findings in many fields. This gel shows a strain of *Salmonella* identified in a food safety experiment.

Specifications

Ordering Information

· · · · · · · · · · · · · · · · · · ·				
Automation Capabilities		Catalog #	Description	
Workflow automated selection	Application driven; user selected or recalled by a protocol	170-8195 Molecular Imager [®] Gel Doc [™] XR+ System with Image Lab Software, PC or Mac, includes darkrd		
Workflow automated execution	Controlled by a protocol via application specific setup for image area, illumination source, filter, analysis, and reporting	Accessories	UV transilluminator, epi-white illumination, camera, cables, Image Lab Software	
Workflow reproducibility	100% repeatability via recallable protocols; from image capture to quantitative analysis and reports	170-8182	XcitaBlue Conversion Screen , includes view goggles; blue conversion screen for viewing SYBR [®] Green, SYBR [®] Safe, GFP, Flamingo, and other	
Autofocus (patent pending)	Precalibrated focus for any zoom setting or sample height	fluorescent gel stains 170-8183 XcitaBlue Conversion Screen and Filter, inc view googles and SVEP® Sofe Filter (170, 9075		
Image flat fielding*	Dynamic; precalibrated and optimized for every application		560DF50); blue conversion screen for viewing SYBR [®] Green, SYBR [®] Safe, and other fluorescent gel stains	
Autoexposure	2 user-defined modes (intense or faint bands)	170-8074	Filter, 520DF30 62 mm, for SYBR® Green/GFP/SYBR® Gold/fluorescein	
Hardware Specifications		170-8075	Filter, 560DF50 62 mm, for SYBR® Safe/Cy3/	
Maximum sample size (L x W)	28 x 36 cm	170-8076	Filter. 630BP30 62 mm. for SYPRO Ruby/Texas Red	
Maximum image area (L x W)	19.4 x 26 cm	170-8077	Filter, 480BP70 62 mm, for Hoechst/coumarin	
Excitation source	Trans-UV and epi-white are standard (302 nm included, with 254 and 365 nm available as options); optional trans-white, self-powered or conversion screen; optional XcitaBlue**	170-6887 170-8097 170-8089 170-7581 170-8184	365 nm UV Lamps, pkg of 6 Standard 302 nm UV Lamps, pkg of 6 Mitsubishi Thermal Printer Mitsubishi Thermal Printer Paper, 4 rolls Gel Alignment Templates, pkg of 3	
Illumination control	3 modes (trans-UV, trans-white, epi-white)	Software 170-9690*	Image Lab Software, PC or Mac, for automated image capture, optimization, and 1-D data analysis	
Detector	CCD	* Included with the imaging system.		
Image resolution	4 megapixels	Acrobat Adaba	and Distances are trademarke of Adaba Custama	
Pixel size (H x V)	4.65 x 4.65 μm	Acrobat, Adobe, and Photoshop are trademarks of Adobe Systems Incorporated. Alexa Fluor, Coomassie Fluor Orange, Qdot, SYBR [®] , SYPRO, and Texas Red are trademarks of Life Technologies Corporation. Cy is a trademark of GE Healthcare group companies. DyLight and Krypton are trademarks of Thermo Fisher Scientific. Exce Microsoft, and PowerPoint are trademarks of Microsoft Corporation. GelGreen and GelRed are trademarks of Biotium, Inc. Hoechst is a trademark of Hoechst GmbH. Mac is a trademark of Apple Inc. Mitsubishi is a trademark of Mitsubishi Companies. Precision Plus Protein Standards are sold under license from Life Technologies Corporation, Carlsbad, CA, for use only by the buyer of the product. The buyer is not authorized to sell or resell this product or its components.		
Filter holder	3 positions (2 for filters, 1 without filter)			
Emission filters	1 included (standard), 3 optional			
Dynamic range	>3.0 orders of magnitude			
Pixel density (gray levels)	4,096			
Instrument size (L x W x H)	36 x 60 x 96 cm			
Instrument weight	32 kg			
Operating Ranges Operating voltage	110/115/230 V AC nominal			
Operating temperature	10–28°C (21°C recommended)			
Operating humidity	<70% noncondensing	Bio-Rad Laboratories, Inc. is licensed by Life Technologies Corporatio		
* U.S. patent 5,951,838.		5,616,502.	oroquera for research use only under 0.5. Faterit NUMDER	

** Use the optional XcitaBlue Kit if performing SYBR® Safe DNA applications; the UV to blue conversion screen allows visualization of DNA samples while protecting user and sample from UV

atent Number SYBR is a trademark of Life Technologies Corporation. Bio-Rad

Laboratories, Inc. is licensed by Life Technologies Corporation to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.



radiation damage.

Bio-Rad Laboratories, Inc.

Life Science Group

Web site www.bio-rad.com USA 800 424 6723 Australia 61 2 9914 2800 Austria 43 1 877 89 01 Belgium 03 710 53 00 Brazil 55 11 3065 7550