

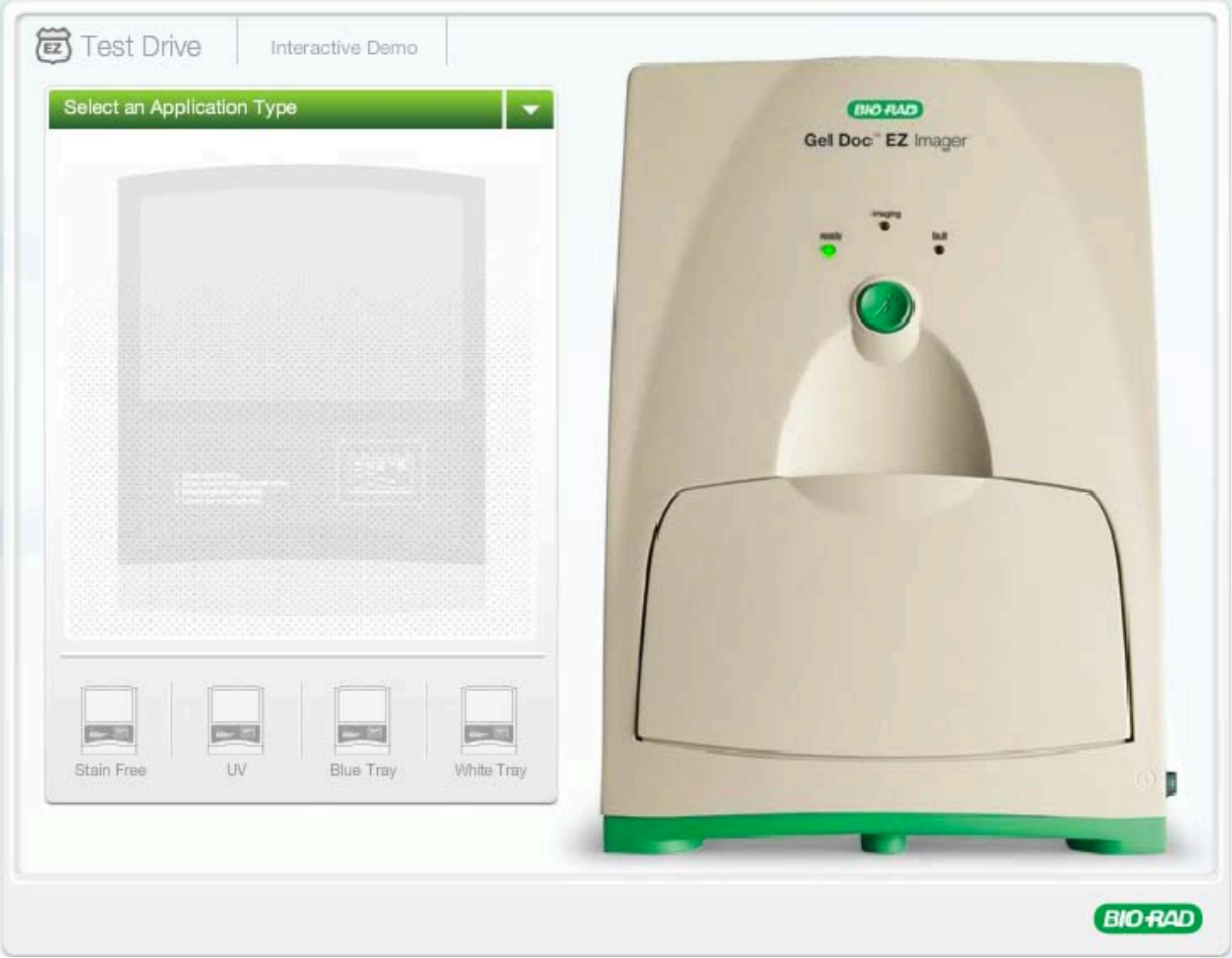


Gel Doc™ EZ Imager

High-quality images and analysis at the touch of a button.

BIO-RAD

Product Demo



Incredibly Smart

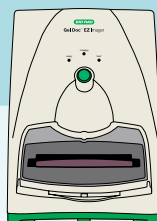
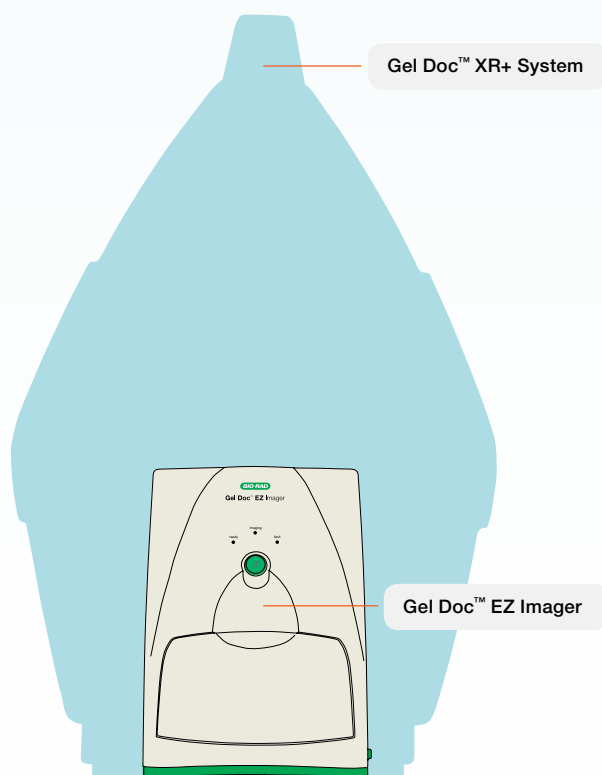
No training required! Just push the green button to get expert images and analysis.

- **Easy to use** — no need for manual control of filters, lens, or lights. Researchers not familiar with the imager can use the system easily
- **High-quality images** — get excellent quality images without any manipulation or user-introduced errors
- **Fast time to results** — eliminate various steps in image acquisition and get the results you need quickly
- **Customizable** — no need to program each and every time. Each tray is customized according to the needs of the individual user

Astoundingly Compact

This imaging system is so small that it can even fit on top of your notebook!

- **Space saving** — free up benchspace, save real estate for running your experiments
- **Convenient** — no need for a dedicated instrument room for your imaging system. Now you can have your imager right next to your electrophoresis apparatus

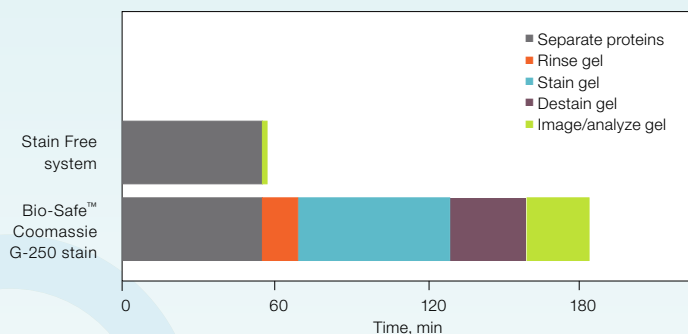
1
Load**2**
Push**3**
Results

38 cm

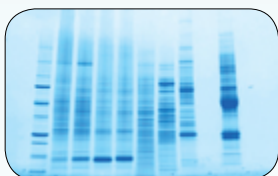
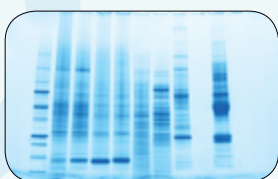


26 cm

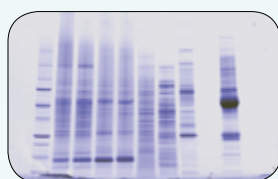
■ Gel Doc EZ System vs. Traditional Coomassie Staining Methods



Stain Free system



Coomassie staining



Expedited Workflow and Increased Free Time Using Stain-Free Technology

Eliminate the traditional SDS-PAGE staining bottleneck using the stain-free method.

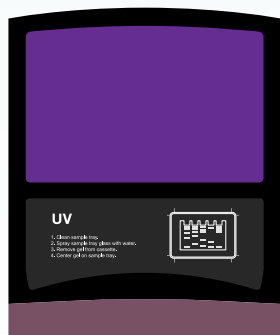
- **Condensed protocol** — convert your 2 hour Coomassie protocol into a 5 minute stain and image step with this stain-free imaging system
- **Compatibility** — Stain-Free gels are western blot compatible, allowing you to check electrophoresis results and quality prior to western blotting
- **Sensitivity** — equal or better sensitivity compared to Coomassie staining
- **Green** — no organic waste disposal concerns

For more information and data, visit www.bio-rad.com/geldocEZ.

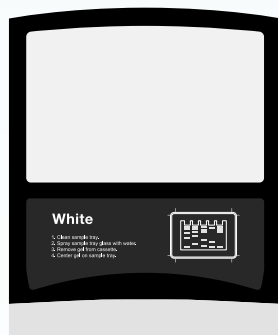
Remarkable Flexibility

Now you have a system that is totally customizable to your needs.

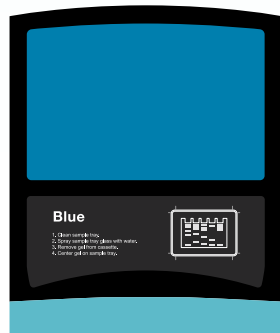
- **Modular design** — use specific trays for specific applications. Clearly defined and color-coded trays eliminate any confusion in usage
- **Flexible options** — purchase only what you want and upgrade when your needs change
- **Simplicity** — create a default protocol once, then use the green button on the front of the instrument to reproducibly use these settings time and again



UV Tray



White Tray



Blue Tray



Stain-Free Tray



Sophisticated Software

The Gel Doc EZ imaging system comes with Image Lab™ software version 3.0, with auto image capture, auto analysis, user preferences, and myriad other features.

- **Completely analyzed results** — no adjusting images, no manual transfer and analysis of data, no plotting, no guesses! Get high-quality images and analyzed results, including relative molecular weights, quantitation of bands, Excel reports, PDFs, and more, within a matter of minutes
- **Reproducibility** — no chances for user-introduced errors. Rely on the system to give consistent results time after time

Picture-Perfect Images Every Time

Obtain high-quality images of your gels every time you push the button.

- **Publication-quality images** — get clean and smooth images that are visually appealing and publication ready
- **Increased image resolution** — decreases pixelation when images are cropped or zoomed
- **Greater functionality** — no need to export images to another program such as Photoshop to change the dpi before importing for publication. You can now define your desired resolution with Image Lab software

For more detailed information and additional sample images, visit www.bio-rad.com/geldocEZ.



Ordering Information

Catalog #	Description
170-8270	Gel Doc EZ System
170-8271	UV Sample Tray
170-8272	White Sample Tray
170-8273	Blue Sample Tray
170-8274	Stain-Free Sample Tray
170-8276	Sample Tray Holder

Coomassie is a trademark of BASF Aktiengesellschaft.

Excel is a trademark of Microsoft Corporation.

Photoshop is a trademark of Adobe Systems Incorporated.

SYBR is a trademark of Invitrogen Corporation.

Bio-Rad Laboratories, Inc. is licensed by Invitrogen Corporation to sell SYPRO products for research use only under U.S. Patent Number 5,616,502.

Krypton is a trademark of Thermo Fisher Scientific Inc.

GelGreen and GelRed are trademarks of Biotium, Inc.

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 20 8732 2339 **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 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500
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 3188 **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 **United Kingdom** 020 8328 2000