

GelDoc Go Imaging System

Go Further





Simple Benchtop Imaging



Select from optimized preset settings for your gels



Click **Acquire** to capture an image



Print or **export** your image in your choice of multiple file types

Ready, Set, GelDoc Go

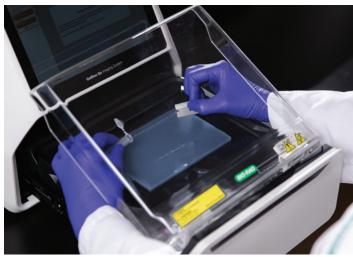
Versatile Applications



Large imaging area in a compact package

Don't let the small footprint fool you — the GelDoc Go Imaging System is compatible with small- and largeformat DNA gels.

For high-throughput applications, capture images of up to four mini gels at once.



Safe gel excision, no matter your stain

Perform gel excision directly on the instrument. The optional UV Shield protects the user when working with traditional DNA stains.

Use the optional Blue Tray for easy direct excision when working with safe stains.

Supported Stains

Protein gels

- Stain-Free gels
- Oriole Fluorescent Gel Stain
- Silver stains
- Coomassie Blue stains
- Coomassie Fluor Orange Stain
- SYPRO Ruby Stain

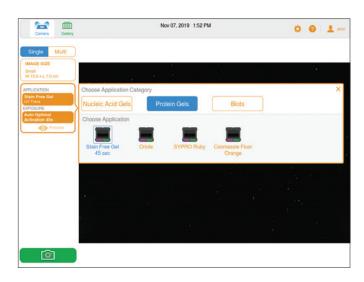
Nucleic acid gels

- Ethidium bromide
- SYBR® Green
- SYBR® Safe
- SYBR® Gold
- GelRed Stain
- GelGreen Stain
- Fast Blast DNA Stain
- UView Stain

Blots

- Stain-Free blots
- Colorimetric blots
- Ponceau S stain

Intelligent Software Packages Designed to Work Together

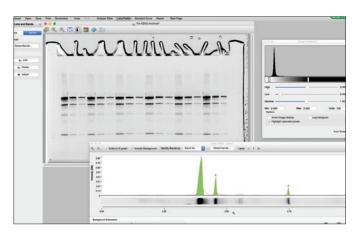


Intuitive image capture with Image Lab Touch Software

Quickly set up the GelDoc Go System to capture your image, add notes, and export or print.

Closely examine your images by pinching and zooming and by adjusting the transform via the touch-screen controls.

Easily manage user accounts and set permissions and security options.



Powerful analysis with desktop Image Lab Software

Image files can be opened in Image Lab Software for accurate, in-depth analysis of your gel data.

Analyzed data can be exported to a spreadsheet or archived in an electronic laboratory notebook.

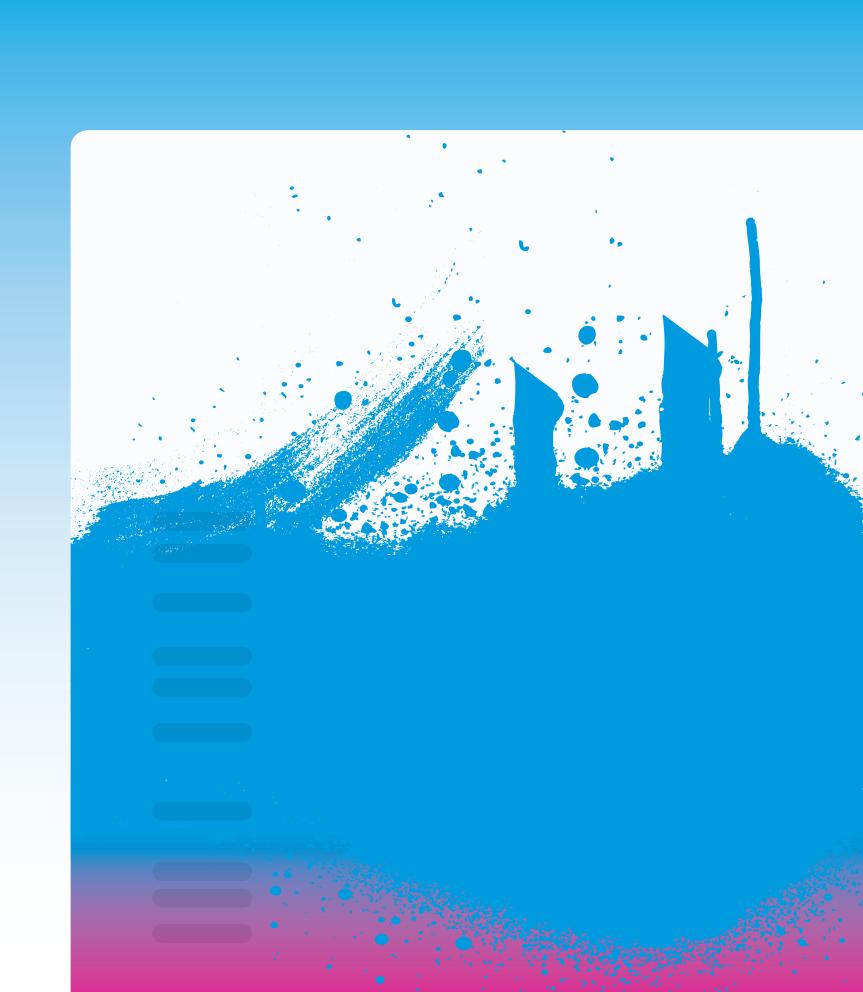
Advanced security options and regulatory compliance features allow you to maintain an unbroken chain of data custody, from capture on the GelDoc Go System through analysis at your computer.

Complete Solutions for Your Needs



Seamless integration into nucleic acid and protein workflows

The GelDoc Go System works with the Bio-Rad products you trust to provide complete solutions for nucleic acid and protein research. Use the C1000 Touch Thermal Cycler and Sub-Cell GT Electrophoresis Cells with the GelDoc Go System to generate and analyze PCR products. For protein workflows, pair the GelDoc Go System with precast gels and Mini-PROTEAN Tetra Vertical Electrophoresis Cells to assess fractions from the NGC Chromatography System.



Specifications

Automation Capabilities

Smart Tray Technology The GelDoc Go Imaging System automatically

recognizes your application-specific tray and adjusts imaging parameters and software options accordingly

Auto-focus Precalibrated focus for any zoom setting Two auto-exposure algorithms (rapid or optimal) Auto-exposure

Hardware Specifications

Touch-screen functionality Multitouch capable

Display resolution 1,024 x 768 pixels

9.7" (24.64 cm) display

21 x 14 cm

Maximum image area

 $(W \times H)$

Detector 6.3 MP CMOS Pixel size 2.4 x 2.4 um Bit depth (gray levels) 65.536

Dynamic range >3.5 orders of magnitude Trans-UVB (standard) Excitation source Epi-white (standard)

> Trans-white (with optional White Tray) Trans-blue (with optional Blue Tray)

Emission filter 535-645 nm (standard)

Data output 16-bit or 8-bit: SCN, TIFF, JPEG image files

Instrument weight ~16 kg (~35 lb) Instrument size (D x W x H) 44.8 x 36.0 x 35.3 cm Operating voltage 100-240 VAC, 50-60 Hz

Operating temperature 10-28°C

Operating humidity 20-80% relative humidity (noncondensing)

Ordering Information

Catalog # Description

Imaging System

12009077 GelDoc Go Imaging System with Image Lab Touch Software

GelDoc Go Imaging System Trays

White Tray for GelDoc Go Imaging System 12012165 12012160 Blue Tray for GelDoc Go Imaging System

12012189 UV/Stain-Free Tray for GelDoc Go Imaging System

Analysis Software

Image Lab Software, on USB drive 12012931 17006130 Image Lab, Security Edition, 1 license 17006171 Image Lab, Security Edition, 5 licenses Image Lab, Security Edition, 10 licenses 17006172

Optional Accessories

UV Shield for GelDoc Go Imaging System 12012164 1708377 Holder for Sample Trays and UV Shield

XcitaBlue Viewing Goggles 1708185 12012190 GelDoc Go Gel Alignment Kit

12012147 GelDoc Go Imaging System IQ/OQ Kit 1708089 Mitsubishi Printer, 100/240 V, USB 1707581 Mitsubishi Thermal Printer Paper, 4 rolls

1703760 Gel Cutter Ruler

Visit bio-rad.com/LetsGo for more information.

BIO-RAD, MINI-PROTEAN, and SUB-CELL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. SYBR is a trademark of Thermo Fisher Scientific Inc. TGX Stain-Free Precast Gels are covered by U.S. Patent Numbers 7,569,130 and 8,007,646. 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 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

Bulletin 7284 Ver A US/EG 19-0703 0420 Sig 0220

