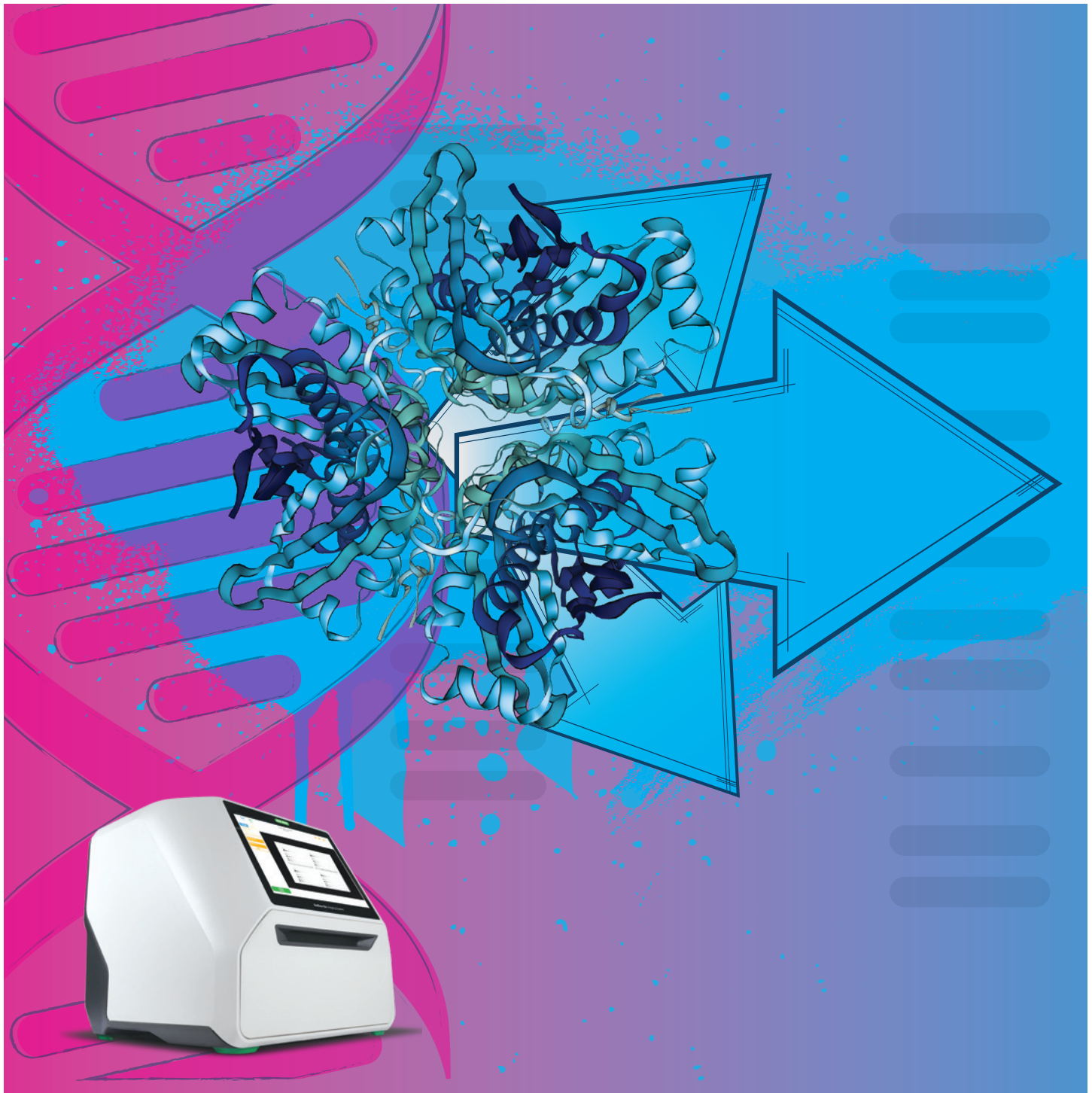


GelDoc Go Imaging System

Go Further



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



Big Impact. Small Footprint.

Benchtop imaging, simplified. All the functionality you expect, with none of the complexity.



White Tray
(Optional)



Blue Tray
(Optional)



UV/Stain-Free Tray
(Included)

Smart Tray Technology automatically adjusts imaging parameters and software options for nucleic acid and protein gels.



Capture sharp and crisp images using Bio-Rad's latest scientific camera technology.

Simple Benchtop Imaging



1 Select from optimized preset settings for your gels



2 Click **Acquire** to capture an image



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

Ready, Set, GelDoc Go

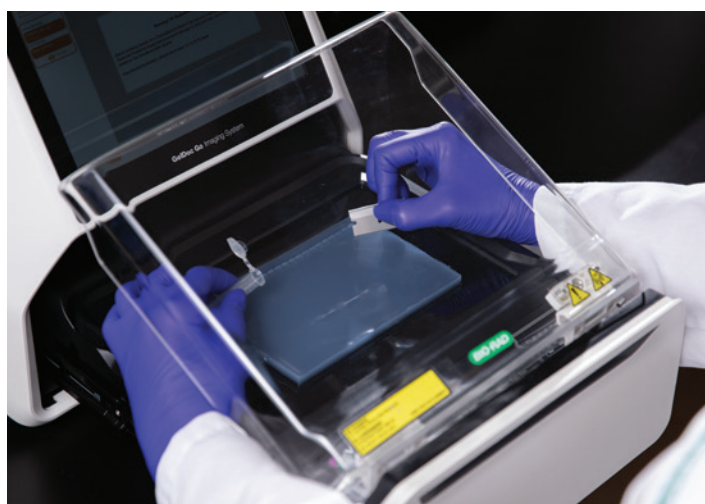
Versatile Applications



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 large-format 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

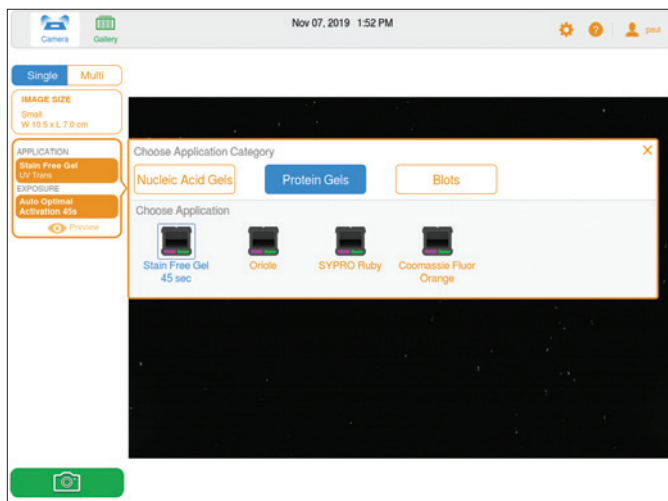
Nucleic acid gels

- 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

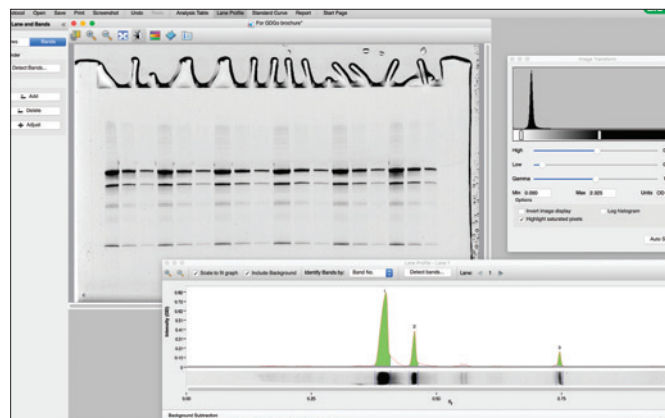


Intuitive image capture with Image Lab Touch Software

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.

A Winning Design

Good Design Australia* has recognized the GelDoc Go System for its exceptional design, stating:

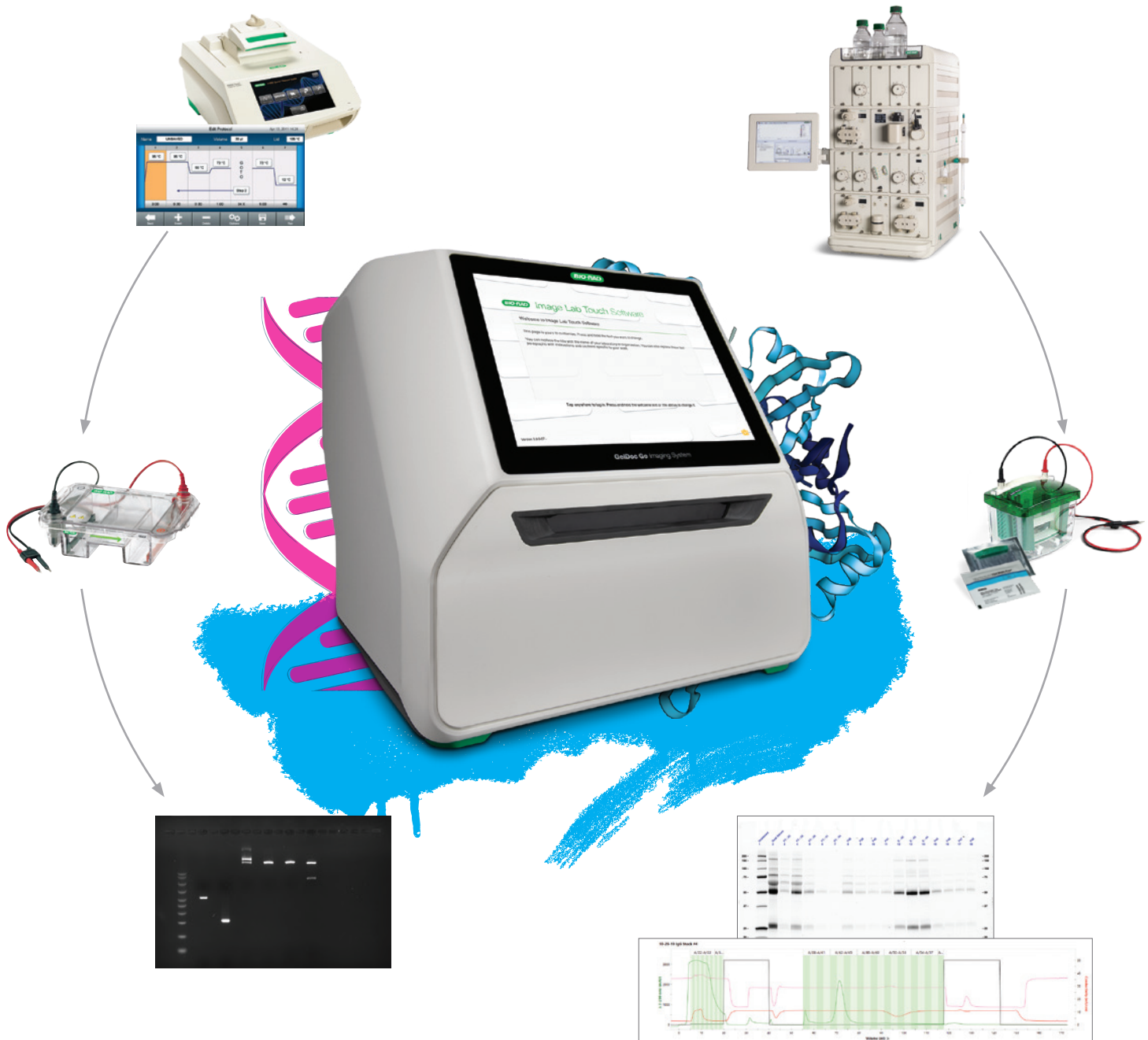


“GelDoc Go is a breakthrough digital imaging instrument used for nucleic acid and protein analysis. Its small footprint, benchtop size, and rapid processing speed make traditional, room-sized instruments redundant, giving medical scientists a powerful tool to improve world health outcomes.”

Design Feature	How It Was Achieved
Streamlined workflow	A telescopic sample drawer provides a single-action opening process, and loading the Smart Tray triggers preset operations and identifies sample types as they are docked
Robust and stable optical chassis	The unique double-skin construction of the die-cast chassis provides stable optical mounting and the bulk of the light sealing, with minimal added weight
Ergonomic	The interface size and angle suit a large range of user heights, giving a clear eyeline and hand access to adjust gels in the cleverly designed forward-leaning drawer front

* Good Design Australia is one of the longest running international design award organizations, promoting excellence in design and innovation since 1958. The Good Design Award, also known as the Good Design Tick, is among the most respected and recognized international design endorsements.

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
Auto-exposure	Two auto-exposure algorithms (rapid or optimal)

Hardware Specifications

Touch-screen functionality	Multitouch capable Display resolution 1,024 x 768 pixels 9.7 in. (24.64 cm) display
Maximum image area (W x H)	21 x 14 cm
Detector	6.3 MP CMOS
Pixel size	2.4 x 2.4 µm
Bit depth (gray levels)	65,536
Dynamic range	>3.5 orders of magnitude
Excitation source	Trans-UVB (standard) 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

12012165	White Tray for GelDoc Go Imaging System
12012160	Blue Tray for GelDoc Go Imaging System
12012189	UV/Stain-Free Tray for GelDoc Go Imaging System

Analysis Software

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

Optional Accessories

12012164	UV Shield for GelDoc Go Imaging System
1708377	Holder for Sample Trays and UV Shield
1708185	XcitaBlue Viewing Goggles
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 Poland 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

