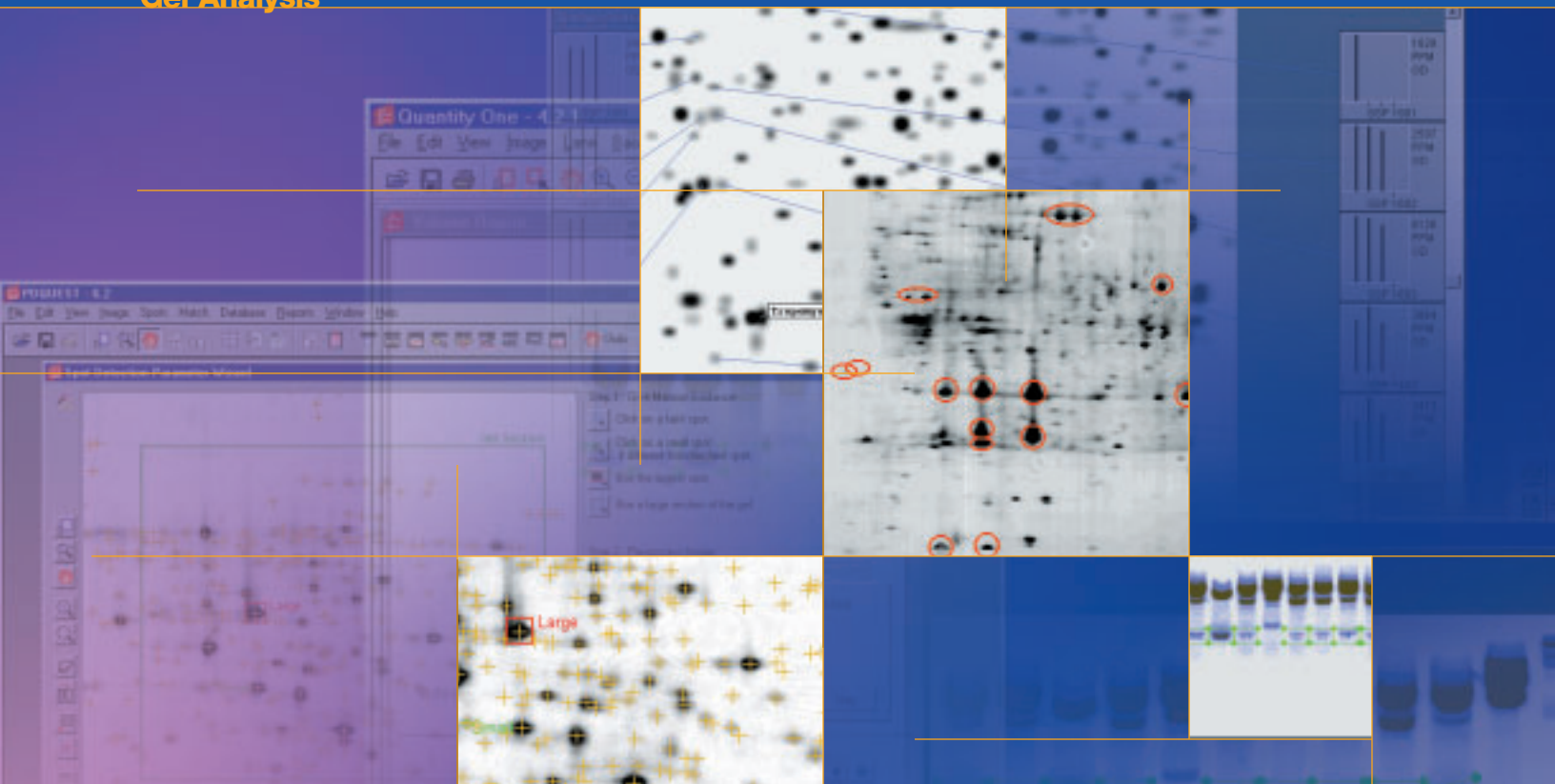


Gel Analysis



The **GS-800**[™] Calibrated Densitometer



Calibrated Quantitation in a Big Way.
Quantitate It All, From Genomics to Proteomics.

BIO-RAD

Calibrated Analysis of Gels, Film, and Blots

Transmissive and Reflective Capabilities

The GS-800 has a built-in calibrated optical density transmissive step tablet for unsurpassed accuracy. Each tablet contains a serial number for complete documentation to an external standard. The ability of the GS-800 to calibrate to an external standard ensures accurate, reproducible results.

29 x 40 cm Format, 12-Bit Precision

If you need to image and quantitate larger gels for enhanced separation of proteins in two dimensions, the GS-800 is the perfect densitometer. It provides 12-bit precision and 36.3 μm resolution to scan and analyze even the closest bands. Unlike conventional desktop scanners, the GS-800 has a large dynamic range, from 0–3.0 OD. This allows accurate quantitation of samples such as Coomassie stained and silver stained gels.

Multiple Wavelengths for Unparalleled Sensitivity

With red, green, and blue color CCD technology, you will easily quantitate most electrophoretic samples or stains. Scan 1-D and 2-D gels, dot and slot blots, films, slides, photographs, and other samples at the proper wavelengths. Moreover, the GS-800 is one of the few densitometers that accommodates wet and thick samples.

Installation Qualification and Operational Qualification (IQ/OQ)

The calibration of the GS-800 can be validated by following the GS-800 Calibrated Densitometer Installation Qualification and Operational Qualification protocols. These protocols are available as an accessory to the GS-800. These straightforward and simple-to-follow IQ/OQ protocols include instructions for installation of the hardware and software, and instructions for verification of the reflectance and transmittance calibration functions.

Windows and Macintosh Software

The GS-800 includes fully integrated Quantity One software for scanning and analysis. For proteomics, Bio-Rad also offers PDQuest 2-D analysis software for 2-D gel matching and database creation (catalog #170-8604 for PC, 170-8612 for Mac). Similarly to Quantity One, PDQuest was designed for full integration with the GS-800.

For more information, contact your local Bio-Rad representative, or visit us on the Web at discover.bio-rad.com.



The ProteomeWorks system is the global alliance between Bio-Rad Laboratories, Inc. and Waters Corporation (Micromass MS Technologies), dedicated to furthering proteomics research.

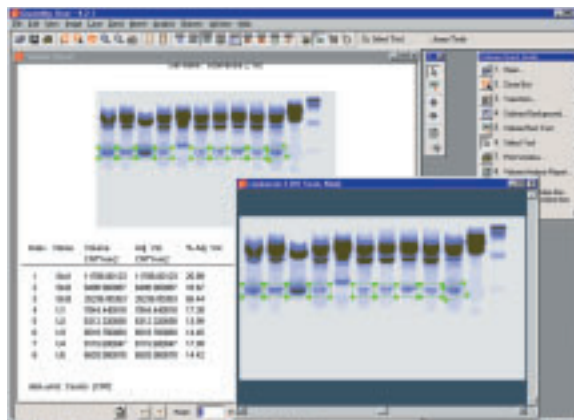


**Bio-Rad
Laboratories, Inc.**

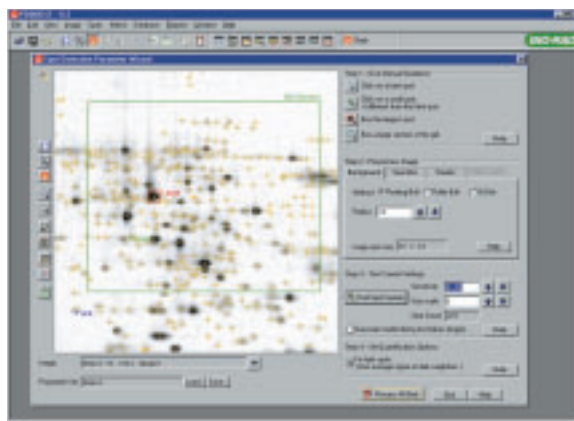
Life Science
Group

Web site www.bio-rad.com **USA** (800) 4BIORAD **Australia** 02 9914 2800 **Austria** (01)-877 89 01 **Belgium** 09-385 55 11 **Brazil** 55 21 507 6191 **Canada** (905) 712-2771 **Czech Republic** + 420 2 41 43 05 32 **China** (86-21) 63052255 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 318 84-177 **Hong Kong** 852-2789-3300 **India** (91-124)-6398112/113/114, 6450092/93 **Israel** 03 951 4127 **Italy** 39 02 216091 **Japan** 03-5811-6270 **Korea** 82-2-3473-4460 **Latin America** 305-894-5950 **Mexico** 52 5 534 2552 to 54 **The Netherlands** 0318-540666 **New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** + 48 22 8126 672 **Portugal** 351-21-472-7700 **Russia** 7 095 721 1404 **Singapore** 65-62729877 **South Africa** 00 27 11 4428508 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717-9555 **Taiwan** (8862) 2578-7189/2578-7241 **United Kingdom** 020 8328 2000

Quantity One® Software



PDQuest™ Software



Specifications

Light source	Fluorescent white
Wavelength	400–750 nm
Operating modes	Transmissive and reflective
Sampling rate	700 dpi
Pixel density	12-bit
Resolution	36.3 μm
Absorbance range	0–3.0 OD
Scanning area	Transmissive: 29 x 40 cm; reflective: 29 x 40 cm
Dimensions	62.6 x 47.3 x 19.3 cm (W x D x H)
Weight	20.8 kg

Ordering Information

Catalog #	Description
170-7980	GS-800 Calibrated Densitometer, PC
170-7981	GS-800 Calibrated Densitometer, Mac
170-7956	GS-800 Calibrated Densitometer Installation Qualification and Operational Qualification
Computers	— please inquire

Coomassie is a trademark of Imperial Chemical Industries PLC. Mac and Macintosh are trademarks of Apple Computer, Inc. Windows is a trademark of Microsoft Corp.