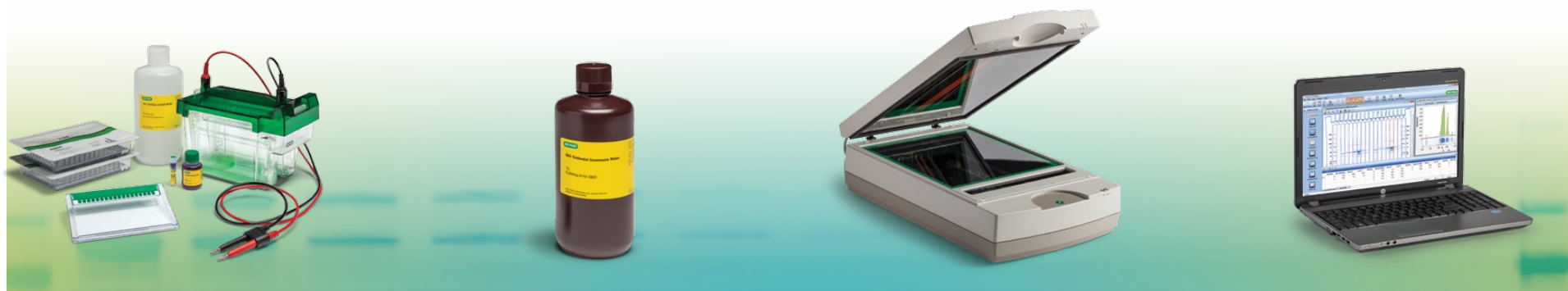


Bio-Rad's Biologics Analysis Workflow™



SEPARATE SAMPLES 1

Faster separations using Criterion™ TGX™ gels

- Gels can be run in as little as 20 minutes
- Load up to 26 samples without decreasing resolution
- Accurate molecular weight estimation with Precision Plus Protein™ standards

Components

- Criterion TGX precast gels
- Criterion™ cell
- Precision Plus Protein standards

STAIN PROTEINS 2

Easier staining with QC colloidal Coomassie stain

- Capable of detecting nanogram levels of protein
- End-point stain ideal for qualitative analysis
- Ready-to-use environmentally friendly stain
- Stain in as little as one hour or as long as overnight

Components

- QC colloidal Coomassie stain

VISUALIZE PROTEINS 3

More reproducible visualization using the GS-900™ calibrated densitometer

- Internal optical density tablet is scanned and used for calibration before each run
- Installation and operation qualification (IQ/OQ) kit confirms the accuracy of the internal target, guaranteeing accurate and reproducible results
- Wide dynamic range permits quantitation of proteins from 0–3.4 OD, ensuring the detection of both highly abundant and dilute proteins

Components

- GS-900 calibrated densitometry system
- IQ/OQ kit

ANALYZE RESULTS 4

Effortless analysis with Image Lab™ software






- Analysis can be automated, saving time and ensuring consistency
- Settings can be tailored to match SOPs
- Tutorials guide users through the analysis process

Components

- Image Lab software
- 21 CFR Part 11 module

BIO-RAD

Ordering Information

Criterion TGX Precast Gels*	 12+2-Well** 45 µl	 18-Well 30 µl	 26-Well 15 µl	 Prep+2-Well 800 µl	 IPG+1-Well 11 cm IPG Strip
7.5% Resolving Gel	567-1023	567-1024	567-1025	—	—
10% Resolving Gel	567-1033	567-1034	567-1035	—	—
12% Resolving Gel	567-1043	567-1044	567-1045	—	—
18% Resolving Gel	567-1073	567-1074	567-1075	567-1072	567-1071
4–15% Linear Gradient	567-1083	567-1084	567-1085	567-1082	567-1081
4–20% Linear Gradient	567-1093	567-1094	567-1095	567-1092	567-1091
8–16% Linear Gradient	567-1103	567-1104	567-1105	567-1102	567-1101
10–20% Linear Gradient	567-1113	567-1114	567-1115	567-1112	567-1111
Any kD™ Gel	567-1123	567-1124	567-1125	567-1122	567-1121

* Criterion TGX gels are sold as single gels. ** Reference well accommodates 15 µl of markers/standards.

Catalog # Description

Protein Standards

- 161-0393 **Precision Plus Protein™ All Blue Standards Value Pack**, 2.5 ml, 250 applications
- 161-0373 **Precision Plus Protein All Blue Standards**, 500 µl, 50 applications
- 161-0396 **Precision Plus Protein™ Unstained Standards Value Pack**, 5 ml, 500 applications
- 161-0363 **Precision Plus Protein Unstained Standards**, 1 ml, 100 applications

Buffers

- 161-0732 **10x Tris/Glycine/SDS**, 1 L
- 161-0747 **4x Laemmli Sample Buffer**, 10 ml
- 161-0737 **2x Laemmli Sample Buffer**, 30 ml
- 161-0710 **2-Mercaptoethanol**, 25 ml

Stain

- 161-0803 **QC Colloidal Coomassie Stain**, 1 L, ready-to-use, nonhazardous colloidal Coomassie G-250 stain for protein polyacrylamide gels

Criterion™ Cell and Systems

- 165-6001 **Criterion Cell**, vertical midi-format electrophoresis cell, includes buffer tank, lid with power cables, 3 sample loading guides (12+2-well, 18-well, 26-well)
- 165-6019 **Criterion Cell and PowerPac™ Basic Power Supply**, 100–120/220–240 V, includes 165-6001 and 164-5050

Instrument

- 170-7991 **GS-900 Calibrated Densitometry System**, PC-compatible calibrated densitometer, cables, Image Lab software, Biologics Analysis Workflow starter kit (Criterion cell, Criterion TGX precast gels, Precision Plus Protein standards, QC colloidal Coomassie stain, buffers)

Regulatory Packages

- 170-7993 **GS-900 Regulatory Tools Package**, GS-900 IQ/OQ kit, set of protocols for installation qualification/operation qualification and NIST-traceable external step tablet for the GS-900 calibrated densitometer, and Image Lab Security Edition, 1 license, 21 CFR Part 11 module
- 170-9692 **Image Lab Security Edition**, 5 licenses, 21 CFR Part 11 module for Image Lab software, use with GS-900, Gel Doc™ EZ, Gel Doc™ XR+, ChemiDoc™ XRS+, and ChemiDoc™ MP imaging systems
- 170-9693 **Image Lab Security Edition**, 10 licenses, 21 CFR Part 11 module for Image Lab software, use with GS-900, Gel Doc EZ, Gel Doc XR+, ChemiDoc XRS+, and ChemiDoc MP imaging systems
- 170-7994 **GS-900 Calibrated Densitometry IQ/OQ Kit**, set of protocols for installation qualification/operation qualification and NIST-traceable external step tablet for the GS-900 calibrated densitometer
- 170-9691 **Image Lab Security Edition**, 1 license, 21 CFR Part 11 module for Image Lab software, use with GS-900, Gel Doc EZ, Gel Doc XR+, ChemiDoc XRS+, and ChemiDoc MP imaging systems



Delivering increased



THROUGHPUT



SENSITIVITY



REPRODUCIBILITY



AUTOMATION

Coomassie is a trademark of BASF Aktiengesellschaft. 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.

BIO-RAD