

STAIN-FREE TECHNOLOGY

Speed to Results with Stain Free

- Fast run times
- Rapid visualization using stain-free enabled imagers
- Long shelf life precast and handcast formats

Utilize Stain-Free Technology in a Variety of Workflows

Fast and Reproducible Protein Separation and Visualization

Bio-Rad's TGX Stain-Free™ gel chemistry uses a unique trihalo compound, which, when activated with UV light, reacts with tryptophan residues allowing for quick visualization of proteins without any staining steps.

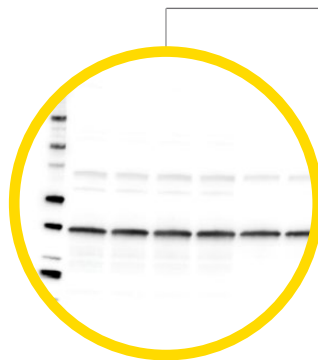
Benefits of Stain-Free Technology

Fast. Run gels in as little as 15 minutes and visualize proteins without staining in 5 minutes.

Low running costs. Suitable for use with inexpensive, standard sample and running buffers.

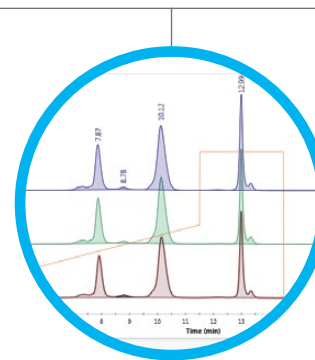
Compatible. Increase speed and efficiency across a variety of workflows.

STAIN-FREE TECHNOLOGY



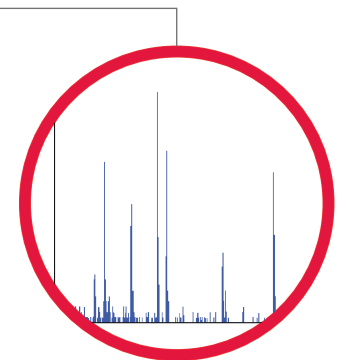
WESTERN BLOTTING

Increase confidence with rapid checkpoints for separation and transfer and get accurate quantitation using total protein normalization.



CHROMATOGRAPHY

Quickly verify protein fraction purity during method development and routine chromatography without any messy, time-consuming staining steps.









MASS SPECTROMETRY

Improve ease of processing and time to results by removing manual staining and destaining steps and allowing for quick visualization of gels.






BIO-RAD

Ordering Information

TGX Stain-Free Precast Gels

Description	 8+1-Well 30 µl	 10-Well 30 µl	 10-Well 50 µl	 12-Well 20 µl	 15-Well 15 µl	 IPG Well 7 cm IPG Strip
Mini-PROTEAN® TGX Stain-Free™ Precast Gels						
7.5% Resolving Gel	456-8029	456-8023	456-8024	456-8025	456-8026	456-8021
10% Resolving Gel	456-8039	456-8033	456-8034	456-8035	456-8036	456-8031
12% Resolving Gel	456-8049	456-8043	456-8044	456-8045	456-8046	456-8041
4–15% Resolving Gel	456-8089	456-8083	456-8084	456-8085	456-8086	456-8081
4–20% Resolving Gel	456-8099	456-8093	456-8094	456-8095	456-8096	456-8091
8–16% Resolving Gel	456-8109	456-8103	456-8104	456-8105	456-8106	456-8101
Any kD™ Resolving Gel	456-8129	456-8123	456-8124	456-8125	456-8126	456-8121

All formats are available in 10-packs (catalog numbers listed) or 2-packs (add an "S" to the end of the catalog number listed).

Description	 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
Criterion™ TGX Stain-Free™ Precast Gels**					
7.5% Gel	567-8023	567-8024	567-8025	—	—
10% Gel	567-8033	567-8034	567-8035	—	—
12% Gel	567-8043	567-8044	567-8045	—	—
18% Gel	567-8073	567-8074	567-8075	567-8072	567-8071
4–15% Gel	567-8083	567-8084	567-8085	567-8082	567-8081
4–20% Gel	567-8093	567-8094	567-8095	567-8092	567-8091
8–16% Linear Gradient	567-8103	567-8104	567-8105	567-8102	567-8101
10–20% Linear Gradient	567-8113	567-8114	567-8115	567-8112	567-8111
Any kD Gel	567-8123	567-8124	567-8125	567-8122	567-8121

* Reference wells accommodate 15 µl of markers/standards. ** Criterion TGX Stain-Free gels are sold as a single gel.

Catalog # Description

TGX Stain-Free™ FastCast™ Acrylamide Solutions and Reagents

161-0181	TGX Stain-Free FastCast Acrylamide Kit, 7.5%
161-0183	TGX Stain-Free FastCast Acrylamide Kit, 10%
161-0185	TGX Stain-Free FastCast Acrylamide Kit, 12%
161-0700	Ammonium Persulfate (APS), 10 g
161-0800	TEMED, 5 ml

Electrophoresis Cells

165-6001	Criterion™ Cell, includes electrophoresis buffer tank, lid with power cables, 3 sample loading guides
165-8004	Mini-PROTEAN® Tetra Cell for Mini Precast Gels, 4-gel vertical electrophoresis system, includes electrode assembly, companion running module, tank, lid with power cables, mini cell buffer dam
165-8000	Mini-PROTEAN Tetra Cell, 4-gel vertical electrophoresis system, 0.75 mm gel thickness
165-8001	Mini-PROTEAN Tetra Cell, 4-gel vertical electrophoresis system, 1.0 mm gel thickness
165-8006	Mini-PROTEAN Tetra Cell, 4-gel vertical electrophoresis system, 1.5 mm gel thickness

Protein Standards

161-0393	Precision Plus Protein™ All Blue Standards Value Pack, 5 pk
161-0396	Precision Plus Protein Unstained Standards Value Pack, 5 pk

Buffers

161-0772	10x Tris/Glycine/SDS, 5 L
161-0737	2x Laemmli Sample Buffer, 30 ml
161-0747	4x Laemmli Sample Buffer, 10 ml

Blotting System

170-4155	Trans-Blot® Turbo™ Starter System, blotting instrument, includes base, 2 cassettes to hold 1–2 midi or up to 4 mini blotting sandwiches, blot roller, and starter consumable kit
----------	--

Detection Reagents

170-5060	Clarity™ Western ECL Substrate, 200 ml
170-5061	Clarity Western ECL Substrate, 500 ml

Catalog # Description

NGC™ Medium-Pressure Chromatography Systems and Modules

788-0001	NGC Quest™ 10 System
788-0003	NGC Quest 10 Plus System, 10 ml/min
788-4004	NGC Sample Pump 100 Module, 100 ml/min
788-4006	NGC Inlet Valve Module
788-4012	NGC Column Switching Valve Module, 10 ml
788-4011	NGC pH Valve Module
788-4010	NGC Buffer Blending Valve Module
788-5017	NGC Air Sensor Module
788-5063	NGC Scout™ Upgrade Kit
788-5064	NGC Discover™ Upgrade Kit

ENrich™ High-Resolution Ion Exchange Columns

780-0001	ENrich Q, 1 ml
780-0021	ENrich S, 1 ml

ENrich High-Resolution Size Exclusion Columns

780-1070	ENrich 70, 25 ml
780-1650	ENrich 650, 25 ml

Imaging Systems

170-8270	Gel Doc™ EZ System with Image Lab™ Software, PC or Mac, includes darkroom, camera, cables, Image Lab software, and stain-free tray. Additional application trays available separately
170-8195	Gel Doc XR+ System with Image Lab Software, PC or Mac, includes darkroom, UV transilluminator, epi-white illumination standard filter, camera, cables, Image Lab software
170-8265	ChemiDoc™ XRS+ System with Image Lab Software, PC or Mac, includes darkroom, UV transilluminator, epi-white illumination standard filter, camera, power supply, cables, Image Lab software
170-8280	ChemiDoc™ MP System with Image Lab Software, PC or Mac, includes darkroom, UV transilluminator, epi-white illumination, standard filter, camera, cables, Image Lab software

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.

Mac is a trademark of Apple, Inc.



**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 11 3065 7550
Canada 905 364 3435 China 86 21 6169 8500 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 9532 220 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300
Israel 03 963 6050 Italy 39 02 216091 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666
New Zealand 64 9 415 2280 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 026 674 55 05
Taiwan 886 2 2578 7189 Thailand 1800 88 22 88 United Kingdom 020 8328 2000