



PROVEN METHOD FOR CONFIRMING PROTEIN PURITY AND YIELD

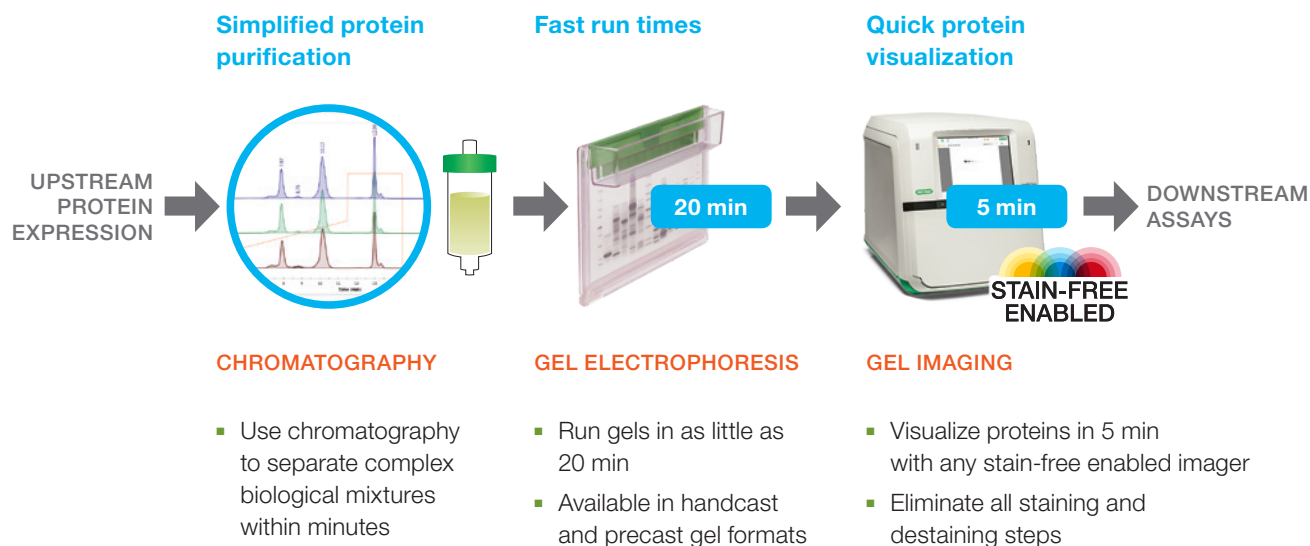
Without the Need for Gel Staining Steps



BIO-RAD

Following chromatography, quickly confirm the purity and yield of your protein with stain-free gels

Stain-free gels provide visual confirmation of protein purity and yield by gel electrophoresis in less than 30 minutes. Stain-free technology is perfect for drug discovery, drug development, and method development for preclinical research.

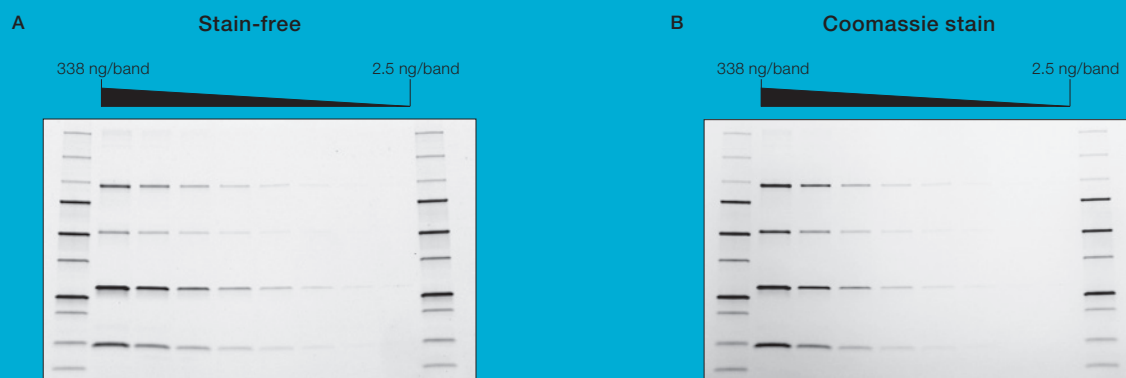


How does stain-free technology work?

Stain-free gel chemistry uses a unique compound, which, when activated, reacts with tryptophan residues in your protein sample to emit a fluorescence signal. This allows the quick visualization of proteins without any staining steps.

Learn more!
Bulletin 6008

Stain-free imaging is as sensitive as Coomassie staining

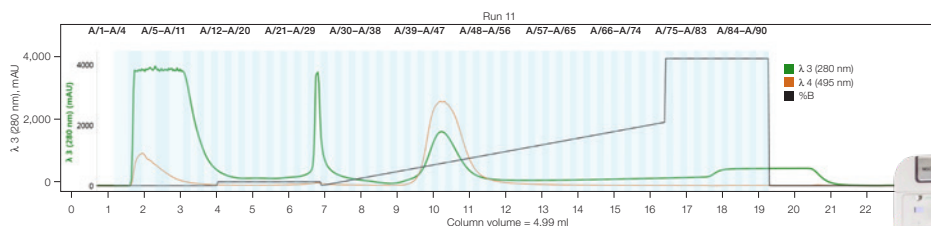


Stain-free and Coomassie-stained gels demonstrate comparable sensitivity when visualizing proteins with varying tryptophan content. 4–20% Criterion TGX Stain-Free Gels were loaded with Precision Plus Protein Unstained Standards and a serial dilution of a protein mixture of four different proteins with varying tryptophan (Trp) content: phosphorylase B (1.4% Trp, 97 kD), glutamic dehydrogenase (0.8% Trp, 56 kD), carbonic anhydrase (2.3% Trp, 31 kD), and lysozyme (3.4% Trp, 14 kD). **A**, gels were run at 300 V for 20 min and immediately imaged with the Gel Doc EZ System; **B**, gels were then stained with Bio-Safe Coomassie Stain and imaged with the Gel Doc EZ System.

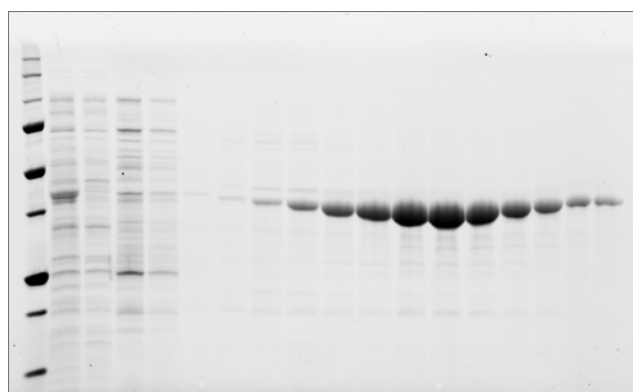
Confidently determine the next steps in your chromatography workflow

Use stain-free gels to quickly:

- Assess protein purity and yield
- Identify which fractions to pool
- Optimize chromatography runs



Learn more!
Bulletin 6530



Visual confirmation of chromatography results in a single stain-free gel.

An 18-well Any kD Criterion TGX Stain-Free Gel was run at 300 V for 20 min. Lane 1 was loaded with 10 μ l Precision Plus Protein Unstained Standards. Lanes 2 and 3 contained 20 μ l of a 1:20 dilution of load and flowthrough, respectively. Lanes 4–18 contained 20 μ l of each fraction.



Stain-free gels are available in both handcast and precast gel formats

FastCast Handcast Gels



- Ready to run in as little as 30 min
- Remain fresh for up to 4 weeks at 4°C
- All the benefits of TGX Stain-Free technology, including fast run times and stain-free visualization of proteins in 5 min

Precast gels



- Available in mini and midi precast gel sizes
- Run up to 26 samples on a single gel
- All the benefits of TGX Stain-Free technology, including fast run times and stain-free visualization of proteins in 5 min



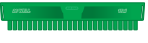


Ordering Information

Mini-PROTEAN TGX Stain-Free Precast Gels

	 10-Well 30 µl	 10-Well 50 µl	 12-Well 20 µl	 15-Well 15 µl	 IPG Well 7 cm IPG Strip
Description					
7.5% Resolving Gel	4568023	4568024	4568025	4568026	4568021
10% Resolving Gel	4568033	4568034	4568035	4568036	4568031
12% Resolving Gel	4568043	4568044	4568045	4568046	4568041
4–15% Resolving Gel	4568083	4568084	4568085	4568086	4568081
4–20% Resolving Gel	4568093	4568094	4568095	4568096	4568091
8–16% Resolving Gel	4568103	4568104	4568105	4568106	4568101
Any kD Resolving Gel	4568123	4568124	4568125	4568126	4568121

All formats are available in 10-packs.

Criterion TGX Stain-Free Precast Gels*

	 12+2-Well** 45 µl	 18-Well 30 µl	 26-Well 15 µl	 Prep+2-Well** 700 µl	 IPG+1-Well** 11 cm IPG Strip
Description					
7.5% Gel	5678023	5678024	5678025	—	—
10% Gel	5678033	5678034	5678035	—	—
12% Gel	5678043	5678044	5678045	—	—
4–15% Gel	5678083	5678084	5678085	5678082	5678081
4–20% Gel	5678093	5678094	5678095	5678092	5678091
8–16% Linear Gradient	5678103	5678104	5678105	5678102	5678101
Any kD Gel	5678123	5678124	5678125	5678122	5678121

* Criterion TGX Stain-Free Gels are sold singly.

** Reference wells accommodate 15 µl of markers/standards.

Catalog # Description

Gel Casting Solutions

1610180	TGX Stain-Free FastCast Acrylamide Starter Kit , 7.5%
1610181	TGX Stain-Free FastCast Acrylamide Kit , 7.5%
1610182	TGX Stain-Free FastCast Acrylamide Starter Kit , 10%
1610183	TGX Stain-Free FastCast Acrylamide Kit , 10%
1610184	TGX Stain-Free FastCast Acrylamide Starter Kit , 12%
1610185	TGX Stain-Free FastCast Acrylamide Kit , 12%

Catalysts

1610700	Ammonium Persulfate (APS) , 10 g
1610800	TEMED , 50 ml

Buffers

1610772	10x Tris/Glycine/SDS , 5 L
1610737	2x Laemmli Sample Buffer , 30 ml
1610747	4x Laemmli Sample Buffer , 10 ml

Electrophoresis Cells

1656001	Criterion Cell , includes electrophoresis buffer tank, lid with power cables, 3 sample loading guides (12+1-well, 18-well, 26-well)
1658004	Mini-PROTEAN Tetra Cell for Mini Precast Gels , 4-gel system includes electrode assembly, mini cell buffer dam, companion running module, tank, lid with power cables
1658000	Mini-PROTEAN Tetra Cell , 10-well, 0.75 mm thickness; 4-gel system includes 5 combs, 1 set of glass plates, 2 casting stands, 4 casting frames, sample loading guide, electrode assembly, mini cell buffer dam, companion running module, tank, lid with power cables
1658001	Mini-PROTEAN Tetra Cell , 10-well, 1.0 mm thickness; 4-gel system includes 5 combs, 1 set of glass plates, 2 casting stands, 4 casting frames, sample loading guide, electrode assembly, mini cell buffer dam, companion running module, tank, lid with power cables
1658006	Mini-PROTEAN Tetra Cell , 10-well, 1.5 mm thickness; 4-gel system includes 5 combs, 1 set of glass plates, 2 casting stands, 4 casting frames, sample loading guide, electrode assembly, mini cell buffer dam, companion running module, tank, lid with power cables

Stain-Free Enabled Imagers

12003153	ChemiDoc Imaging System , blot and gel imaging system, UV/visible light imaging, chemiluminescence, upgradeable for multiplex fluorescence detection; includes internal computer, 12" touch-screen display, Image Lab Touch Software, blot/UV/stain-free sample tray
12003154	ChemiDoc MP Imaging System , blot and gel imaging system, UV/visible light imaging, chemiluminescence, 5 fluorescence channels (red, green, and blue [RGB], far red, near infrared); includes internal computer, 12" touch-screen display, Image Lab Touch Software, blot/UV/stain-free sample tray
1708265	ChemiDoc XRS+ System with Image Lab Software , PC or Mac, includes darkroom, UV transilluminator, epi-white illumination standard filter, camera, power supply, cables
1708270	Gel Doc EZ System , PC or Mac, includes Image Lab Software, darkroom, camera, cables
1708195	Molecular Imager Gel Doc XR+ System with Image Lab Software , PC or Mac, includes darkroom, UV transilluminator, epi-white illumination standard filter, camera, cables

Mini-PROTEAN and Criterion Empty Cassettes and Combs

4560001	Mini-PROTEAN Empty Cassettes , 1.0 mm thick, IPG well, pkg of 50
4560003	Mini-PROTEAN Empty Cassettes , 1.0 mm thick, 10-well, pkg of 50
4560005	Mini-PROTEAN Empty Cassettes , 1.0 mm thick, 12-well, pkg of 50

Catalog # Description

4560006	Mini-PROTEAN Empty Cassettes , 1.0 mm thick, 15-well, pkg of 50
4560011	Mini-PROTEAN Combs , IPG well, pkg of 50, for use with 4560001
4560013	Mini-PROTEAN Combs , 10-well, pkg of 50, for use with 4560003
4560015	Mini-PROTEAN Combs , 12-well, pkg of 50, for use with 4560005
4560016	Mini-PROTEAN Combs , 15-well, pkg of 50, for use with 4560006
1653310	Spacer Plates with 0.75 mm Integrated Spacers , 5
1653311	Spacer Plates with 1.0 mm Integrated Spacers , 5
1653312	Spacer Plates with 1.5 mm Integrated Spacers , 5
3459901	Criterion Empty Cassettes , 1 mm, 12+2-well, pkg of 10
3459902	Criterion Empty Cassettes , 1 mm, 18-well, pkg of 10
3459903	Criterion Empty Cassettes , 1 mm, 26-well, pkg of 10
3459904	Criterion Empty Cassettes , 1 mm, prep+2-well, pkg of 10
3459906	Criterion Empty Cassettes , 1 mm, IPG+1-well, pkg of 10
Protein Standards	
1610393	Precision Plus Protein All Blue Prestained Standards Value Pack , 5 x 500 µl, 250 applications
1610396	Precision Plus Protein Unstained Standards Value Pack , 5 x 1,000 µl, 500 applications
NGC Medium-Pressure Chromatography Systems and Modules	
7880001	NGC Quest 10 System
7880003	NGC Quest 10 Plus System
7880002	NGC Quest 100 System
7880004	NGC Quest 100 Plus System
7880005	NGC Scout 10 System
7880007	NGC Scout 10 Plus System
7880009	NGC Discover 10 System
7884010	NGC Buffer Blending Valve Module
7884006	NGC Inlet Valve Module
7884011	NGC pH Valve Module
7884012	NGC Column Switching Valve Module , 10 ml
7884026	NGC Column Switching Valve Module , 100 ml
7884013	NGC Outlet Valve Module
7410002	BioFrac Fraction Collector
17002070	NGC Fraction Collector
ENrich High-Resolution Ion Exchange Columns	
7800001	ENrich Q , 1 ml
7800003	ENrich Q , 8 ml
7800021	ENrich S , 1 ml
7800023	ENrich S , 8 ml
ENrich High-Resolution Size Exclusion Columns	
7801070	ENrich 70 , 24 ml
7801650	ENrich 650 , 24 ml

Bio-Rad, Mini-PROTEAN, and Molecular Imager are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. TGX Stain-Free Precast Gels are covered by U.S. Patent Numbers 7,569,130 and 8,007,646. 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.

All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00
Germany 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600
Japan 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30
Poland 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580
Sweden 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300
United Kingdom 44 01923 47 1301

