

Comparison of Prestained Protein Standards

Bio-Rad's Precision Plus Protein™ Prestained Standards Offer the Sharpest Results



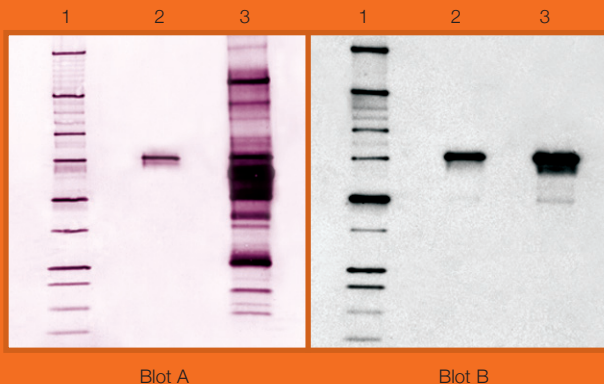
Prestained standards were run on a Bio-Rad Criterion™ 4–20% precast gel in Tris-glycine-SDS running buffer at 200 V for 45 min. Molecular weights shown are for Precision Plus Protein standards.

- 1 Bio-Rad** Precision Plus Protein dual color standards
- Invitrogen BenchMark prestained protein ladder
- Invitrogen SeeBlue Plus2 prestained protein standard
- Amersham Biosciences Rainbow full range molecular weight markers
- 5 Bio-Rad** Precision Plus Protein dual color standards
- 6 Bio-Rad** Precision Plus Protein dual color standards
- Sigma high range color markers
- BMA ProSieve color protein markers
- Invitrogen MultiMark multicolored standard
- 10 Bio-Rad** Precision Plus Protein dual color standards

BIO-RAD

Precision Plus Protein Unstained Standards

Precise Molecular Weight Determinations on Western Blots



Precision Plus Protein standards produce clean signals with sensitive detection by chemiluminescence.

Duplicate Criterion 4–20% Tris-glycine gels were transferred onto nitrocellulose using the Criterion blotter for 30 min at 100 V. Lane 1, Precision Plus Protein unstained standards; lane 2, transferrin; lane 3, total human serum.

Blot A was stained with colloidal gold total protein stain. Blot B was incubated with anti-transferrin followed by StrepTactin-HRP for detection of the Precision Plus Protein standards and GAR-HRP for detection of transferrin. The bands were visualized with Bio-Rad's Immuno-Star™ HRP chemiluminescent detection kit.

Ordering Information

Catalog #	Description
161-0374	Precision Plus Protein Dual Color Standards, 50 applications
161-0373	Precision Plus Protein All Blue Standards, 50 applications
161-0363	Precision Plus Protein Unstained Standards, 100 applications
161-0380	Precision Protein™ StrepTactin-HRP conjugate, 0.3 ml
161-0382	Precision Protein StrepTactin-AP conjugate, 0.3 ml

StrepTactin is a trademark of Institut für Bioanalytik GmbH.

NOTICE TO PURCHASER: LIMITED LICENSE

This product is 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

**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 31 3689 6600 **Canada** 905 364 3435 **China** 86 20 8732 2339 **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** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500 **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** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000