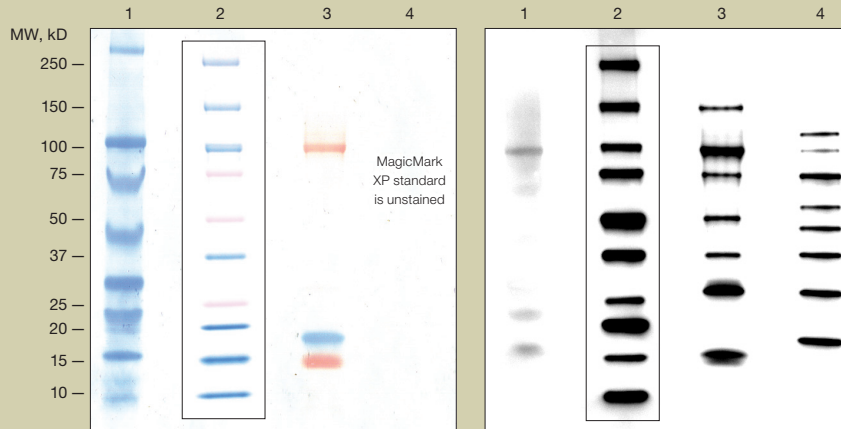
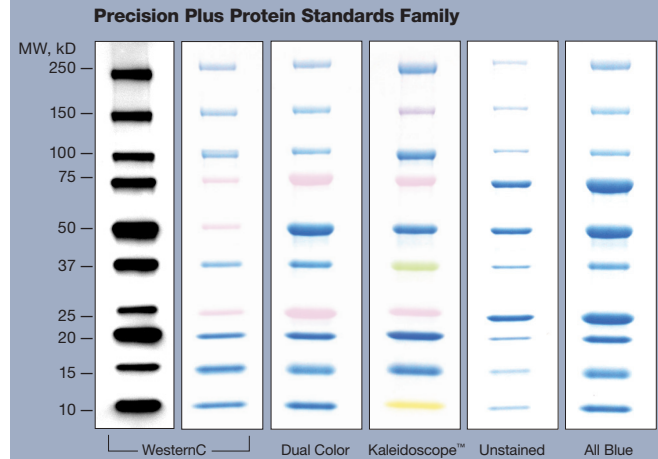


Comparison of Chemiluminescent Protein Standards

Precision Plus Protein™TM WesternC™ standards are prestained chemiluminescent protein standards. The Precision Plus Protein WesternC standards feature the same sharp, clear bands as the other standards in the Precision Plus Protein standards family.



Chemiluminescent standards run on a Criterion™ 4–20% precast gel in Tris/glycine/SDS running buffer at 200 V for 60 min (left) and on a chemiluminescent western blot (right). Molecular weights shown are for Precision Plus Protein WesternC standard bands. Lane 1, Pierce chemiluminescent BlueRanger marker mix; lane 2, Bio-Rad Precision Plus Protein WesternC standards; lane 3, GE Healthcare ECL DualVue western blotting markers; lane 4, Invitrogen MagicMark XP western protein standard.



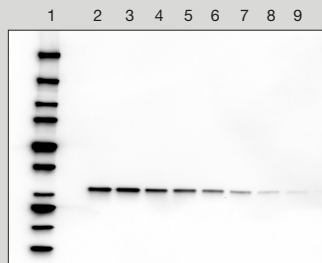
The Precision Plus Protein standards family offers extremely accurate and consistent protein standards for electrophoresis and western blotting experiments. All Precision Plus Protein standards generate a standard curve with R^2 values >0.99 , demonstrating a strong linear relationship between the proteins' log molecular weight and relative migration distance on a gel. The five products in the Precision Plus Protein standards family offer choices and solutions for all your molecular weight estimation needs.

BIO-RAD



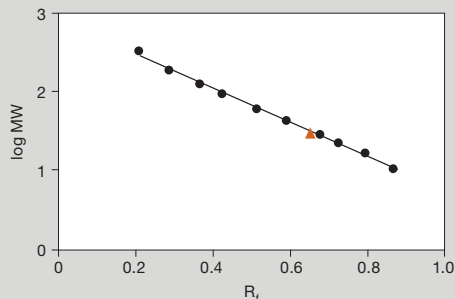
Outstanding Substrate for Charge-Coupled Device (CCD) Imaging

Immun-Star™ WesternC™ chemiluminescent detection kit produces strong, long-lasting signals for CCD imaging.



Western blot detection of 27 kD protein and Precision Plus Protein WesternC standards using the Immun-Star WesternC chemiluminescent detection kit.

Maximum sensitivity achievable with the Immun-Star WesternC kit is in the mid-femtogram range. A gel run with 5 μ l of Precision Plus Protein WesternC standards (lane 1) and a dilution series of *E. coli* lysate containing an overexpressed 27 kD protein (lanes 2–9) was transferred to a nitrocellulose membrane. The dilutions are: 200 ng, 150 ng, 100 ng, 75 ng, 50 ng, 25 ng, 12 ng, 6 ng. The blot was probed with a primary antibody specific for the 27 kD protein, then incubated with StrepTactin-horseradish peroxidase (HRP) and a secondary antibody conjugated to HRP. After incubation in the Immun-Star WesternC detection solution for 5 min, the blot was imaged using the Molecular Imager® ChemiDoc™ XRS system.



Western blot analysis using Bio-Rad Quantity One® 1-D analysis software.

Molecular weight of an unknown protein is estimated using a standard curve generated by Precision Plus Protein WesternC standards. The curve was produced by plotting the log molecular weight (MW) versus the relative migration distance (R_f) of each protein standard through an SDS-PAGE gel. The R^2 value of 0.997 shows the strong linearity of Precision Plus Protein WesternC standards and the accurate estimation of the molecular weight of the 27 kD unknown protein (\blacktriangle). The blot was developed using the Immun-Star WesternC chemiluminescent kit and imaged with the Molecular Imager ChemiDoc XRS system.

Ordering Information

Catalog #	Description
170-5070	Immun-Star WesternC Chemiluminescent Kit , includes 50 ml luminol/enhancer, 50 ml peroxide solution
170-5046	Immun-Star Goat Anti-Rabbit (GAR)-HRP Conjugate , 2 ml, lyophilized
170-5047	Immun-Star Goat Anti-Mouse (GAM)-HRP Conjugate , 2 ml, lyophilized
161-0376	Precision Plus Protein WesternC Standards , 250 μ l, 50 applications
161-0385	Precision Plus Protein WesternC Pack , 50 applications of WesternC standards and 50 applications of StrepTactin HRP
170-8650	Molecular Imager® VersaDoc™ MP 5000 System , PC or Mac, 100/240 V, includes CCD camera, darkroom, power supply, cables, epi-illumination, transillumination, fluorescent reference plate, focusing target, Quantity One software, instructions
170-8070	Molecular Imager ChemiDoc XRS System , PC, includes darkroom, UV transilluminator, epi-white illumination, fluorescent reference plate, camera, power supply, PCI digitizing card, cables, Quantity One software, instructions
170-8071	Molecular Imager ChemiDoc XRS System, Mac
161-0380	Precision Protein™ StrepTactin-HRP Conjugate , 300 μ l, 150 applications
161-0381	Precision Protein StrepTactin-HRP Conjugate , 125 μ l, 50 applications

BlueRanger is a trademark of Pierce Chemical Company. ECL DualVue is a trademark of GE Healthcare. Mac is a trademark of Apple Inc. MagicMark is a trademark of Invitrogen Corporation. StrepTactin is a trademark of Institut für Bioanalytik GmbH. StrepTactin is covered by German patent application P 19641876.3. Bio-Rad Laboratories, Inc. is licensed by Institut für Bioanalytik GmbH to sell these products for research use only.

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.