

Protein Standards Selection Guide

From the Leader in Protein Standards



	Precision Plus Protein™ Standards						Prestained Natural Standards				Unstained Natural Standards				Specialty Standards		
	WesternC™	Kaleidoscope™	Dual Xtra	Dual Color	All Blue	Unstained	Broad Range	Low Range	High Range	Natural Kaleidoscope	Broad Range	Low Range	High Range	Polypeptide	IEF	2-D	Standard Plugs***
Molecular weight (MW)/ pI range	10–250 kD	10–250 kD	2–250 kD	10–250 kD	10–250 kD	10–250 kD	6.9–210 kD	14–97 kD	45–200 kD	7.6–216 kD	6.5–200 kD	14–97 kD	45–200 kD	1.4–26.6 kD	4.45–9.6 pI	17.5–76 kD 4.5–8.5 pI	10–250 kD
Number of proteins	10	10	12	10	10	10	8	6	4	7	9	6	5	6	9	7	10
Electrophoresis																	
SDS-PAGE	•	•	•	•	•	•	•	•	•	•	•	•	•	•	-	-	•
Accurate MW estimation	•	•	•	•	•	•	-	-	-	-	•	•	•	•	-	-	•
Multicolored	•	•	•	•	-	-	-	-	-	•	-	-	-	-	-	-	-
Coomassie staining	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•
Fluorescent staining	-	-	-	-	-	•	-	-	-	-	•	•	•	•	-	•	•
2-D electrophoresis	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	•	-
IEF	-	-	-	-	-	-	-	-	-	-	-	-	-	-	•	-	-
Plug format for use in gels with no reference well	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	-	•
Blotting																	
Monitoring transfer efficiency	•	•	•	•	•	-	•	•	•	•	-	-	-	-	-	-	-
Coomassie staining	•	•	•	•	•	•	•	•	•	•	•	•	•	•	-	•	•
Fluorescent staining	-	-	-	-	-	•	-	-	-	-	•	•	•	•	-	•	•
Fluorescent blotting*	•	•	•	•	•	-	-	-	-	-	-	-	-	-	-	-	-
Immunodetection**	•	-	-	-	-	•	-	-	-	-	-	-	-	-	-	-	•
Catalog Numbers																	
Single unit	161-0385	161-0375	161-0377	161-0374	161-0373	161-0363	161-0318	161-0305	161-0309	161-0324	161-0317	161-0304	161-0303	161-0326	161-0310	161-0320	161-0378
Value pack of 5 units	161-0398	161-0395	161-0397	161-0394	161-0393	161-0396	-	-	-	-	-	-	-	-	-	-	-

* These standards have fluorescent properties and can be used for fluorescent blotting applications. See bulletin 5723 for details on using Precision Plus Protein WesternC standards for fluorescent multiplexing. Precision Plus Protein Dual Xtra standards (161-0377) are recommended for fluorescent blot analysis of proteins between 5 and 250 kD.

** Immunodetection via addition of a Precision Protein™ StrepTactin and horseradish peroxidase (HRP) or StrepTactin and alkaline phosphatase (AP) conjugate, which will bind to the internal Strep-tags on the proteins.

*** 24 unstained plugs for 2-D gels.



Protein Standards for Electrophoresis and Western Blotting

Recombinant Precision Plus Protein Standards

Prestained

Unstained

Natural Standards

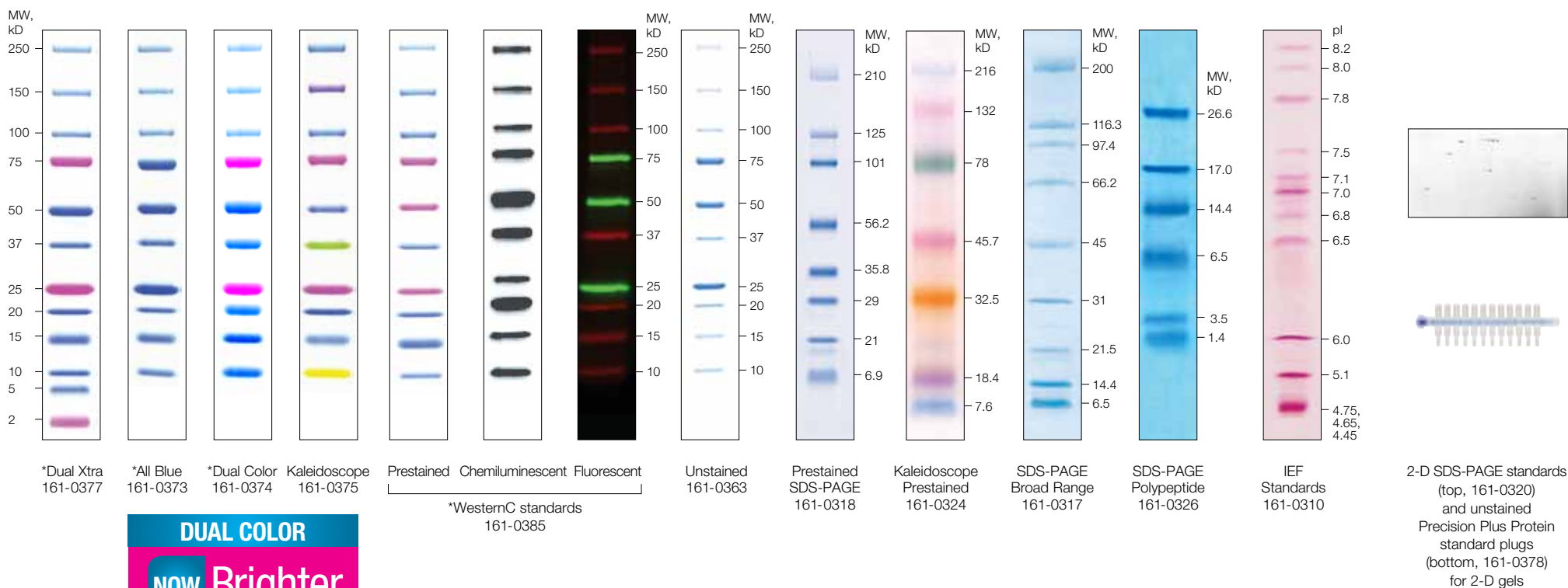
Prestained

Unstained

Specialty Standards

Prestained

Unstained



* These standards have fluorescent properties and can be used for fluorescent blotting applications. See bulletin 5723 for details on using Precision Plus Protein WesternC standards for fluorescent multiplexing.

Note: Value packs of 5 are available for the following Precision Plus Protein standards: All Blue (161-0393), Dual Color (161-0394), Kaleidoscope (161-0395), Dual Xtra (161-0397), WesternC (161-0398), and Unstained (161-0396)

For more information, please visit us on the web at www.bio-rad.com/proteinstandards.

Strep-tag technology for western blot detection is covered by U.S. Patent Number 5,506,121 and by UK Patent Number 2,272,698.

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.

