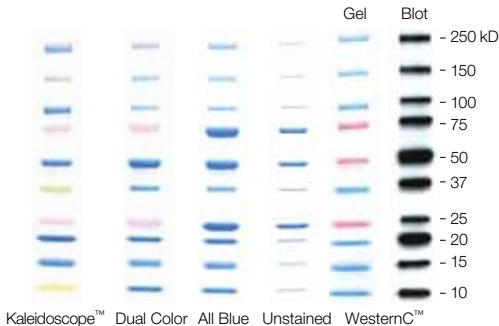


Did You Know?

Bio-Rad offers a complete family of Precision Plus Protein™ standards with the same migration pattern for all your protein electrophoresis and blotting needs.



BIO-RAD

Ordering Information

Catalog #	Description
161-0373	Precision Plus Protein All Blue Standards , 500 μ l, 50 applications
161-0374	Precision Plus Protein Dual Color Standards , 500 μ l, 50 applications
161-0375	Precision Plus Protein Kaleidoscope Standards , 500 μ l, 50 applications
161-0363	Precision Plus Protein Unstained Standards , 1 ml, 100 applications
161-0385	Precision Plus Protein WesternC Pack (standards and conjugate) , 50 applications
161-0376	Precision Plus Protein WesternC Standards , 250 μ l, 50 applications

For more information, go to:

www.bio-rad.com/ppstandards/.

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.