
Introducing Precision Plus Protein™ Standards

Available Now!

The reliability of Precision Protein™ standards, plus two colors and even greater accuracy!

BIO-RAD

The Precision Protein™ standards that you have relied on for superior performance have been made even better!

You know that Precision Protein standards deliver the sharpest bands and molecular weight accuracy. Precision Plus Protein standards deliver the same reliability, plus:

Instant Orientation

Three reference bands in the Precision Plus Protein standards provide instant orientation of gels and blots for easy analysis.

Even Higher Accuracy

Precision Plus Protein standards provide an additional protein band at 20 kD for more accurate estimation of molecular weights in the 10–20 kD range. Also, a new 15 kD protein band that provides a more accurate plot ($R^2 > 0.99$) of migration vs. log molecular weight in this size range has been substituted for the previous protein.

Two-Color Convenience

Precision Plus Protein standards are also offered in a dual color, prestained blue and pink version. Now you can monitor the progress of your gel electrophoresis more easily and quickly focus on the region of interest.

Get the Benefits of These New, Improved Standards the Next Time You Order!

Get even greater accuracy and instant orientation, plus the Precision Protein features you rely on.

True Molecular Weight

Each protein in the unstained standards was designed for ultimate molecular weight accuracy. Molecular weight is confirmed by mass spectrometry.

Clean, Sharp Bands

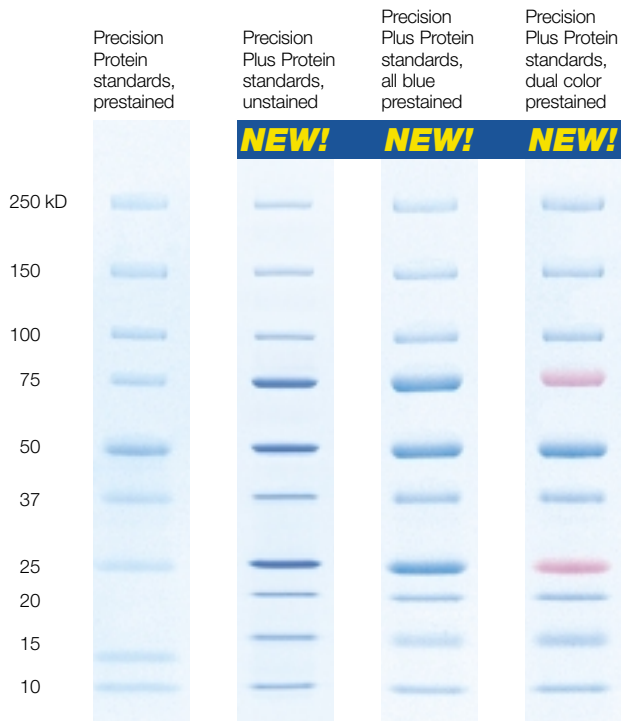
Ten highly purified recombinant proteins are blended to provide the sharpest protein markers available.

Reproducibility

Proprietary staining technology provides batch-to-batch molecular weight consistency and reproducible electrophoretic migration in the prestained standards.

Blot Detection

The *Strep*-tag sequence in the unstained standards enables parallel detection on western blots.



Try Precision Plus Protein Standards—Today!

Ordering Information

Catalog #	Description
161-0363	Precision Plus Protein Standards, unstained, 100 applications per tube
161-0373	Precision Plus Protein Standards, all blue prestained, 50 applications per tube
161-0374	Precision Plus Protein Standards, dual color prestained, 50 applications per tube

If you would like to try these standards and compare them to what you are using now, order a free sample (while supplies last):

161-0363S	Precision Plus Protein Standards, unstained, 50 μ l sample
161-0373S	Precision Plus Protein Standards, all blue prestained, 50 μ l sample
161-0374S	Precision Plus Protein Standards, dual color prestained, 50 μ l sample

The molecular weights of the Precision Plus Protein all blue and dual color prestained standards are confirmed by migration in a Laemmli SDS-PAGE system. For electrophoretic determination of molecular weights at an accuracy >95%, unstained Precision Plus Protein standards (catalog #161-0363) should be used.

Strep-tag is a trademark of Institut für Bioanalytik GmbH. *Strep-tag* technology for western blot detection is covered by US patent 5,506,121 and UK patent 22 72 698. Bio-Rad is licensed to sell these products for research use only.

Visit us on the Web at **www.bio-rad.com**.
Call toll free at 1-800-4BIORAD (1-800-424-6723);
outside the US, contact your local sales office.

The Bio-Rad logo consists of the words "BIO-RAD" in a bold, white, sans-serif font. The text is centered within a dark blue rounded rectangular background.