

PageRuler Prestained Protein Ladder

26616 26617 26618

2353.1

Number	Description
26616	PageRuler Prestained Protein Ladder, 2 × 250μL
26617	PageRuler Prestained Protein Ladder, 10 × 250μL
26618	PageRuler Prestained Protein Ladder, 25μL

Storage Buffer: 62.5mM Tris•H₃PO₄ (pH 7.5 at 25°C), 1mM EDTA, 2% (w/v) SDS, 10mM DTT, 1mM NaN₃ and 33% (v/v) glycerol.

Storage: Upon receipt store at -20°C. Product is shipped with an ice pack.

Introduction

The Thermo Scientific™ PageRuler™ Prestained Protein Ladder is a prestained mixture of ten recombinant proteins ranging from 10K to 180K. Three different chromophores are bound to the proteins, producing a brightly colored ladder (see website for product images). The protein ladder is conveniently packaged and ready to use with no heating, diluting or additional reducing agent necessary.

Important Product Information

- Do not boil the protein ladder.
- Store the protein ladder for up to three months at 4°C or for one year at -20°C.
- The molecular weights of the proteins have a lot-to-lot variation of approximately 5%.
- In low-percentage gels (< 10%), the low-molecular weight proteins in the ladder may migrate with the dye front.
- The large proteins (> 100K) in the ladder may require longer transfer times or higher transfer voltages for Western blotting.
- The mobility of prestained proteins can vary in different SDS-PAGE buffer systems; however, they are suitable for approximate molecular weight determination when calibrated against unstained standards in the same system. See website for migration patterns in different electrophoresis conditions.

Procedure for Use in Polyacrylamide Gel Electrophoresis

1. Thaw the ladder at room temperature. Do not boil protein ladder.
2. Mix gently and thoroughly to ensure that the solution is homogeneous.
3. Load an appropriate volume of the ladder onto the gel.
 - Mini-gel: 5μL per well (0.75-1.0mm thick) or 10μL per well (1.5mm thick)
 - Large gel: 10μL per well (0.75-1.0mm thick) or 20μL per well (1.5mm thick)
4. Return the unused protein ladder to -20°C for up to one year or 4°C for up to three months.

Related Thermo Scientific Products

Please see the catalog or website for a complete listing of protein gels and Western blotting products.

26610	Unstained Protein Molecular Weight Marker, 2 × 1mL
26612	Prestained Protein Molecular Weight Marker, 2 × 250µL
26614	PageRuler™ Unstained Protein Ladder, 2 × 250µL
26619	PageRuler™ Plus Prestained Protein Ladder, 2 × 250µL
26630	PageRuler™ Broad Range Unstained Protein Ladder, 2 × 250µL
26632	PageRuler™ Low Range Unstained Protein Ladder, 2 × 250µL
26634	Spectra™ Multicolor Broad Range Protein Ladder, 2 × 250µL
26625	Spectra™ Multicolor High Range Protein Ladder, 2 × 250µL
26628	Spectra™ Multicolor Low Range Protein Ladder, 250µL
84785	SuperSignal™ Molecular Weight Protein Standards, 250µL
84786	SuperSignal™ Enhanced Molecular Weight Protein Standards, 250µL
25244	Precise™ Protein Gels, 4-20%, 15 well
84713	Pierce™ Protein Gels, 4-20%, 12 well
24615	Imperial™ Protein Stain, 1L
24594	GelCode™ Blue Safe Protein Stain, 1L

General References

- Alegria-Schaffer, A., *et al.* (2009). Performing and optimizing Western blots with an emphasis on chemiluminescent detection. *Methods Enzymol* **463**:573-99.
- Burnette, W.N. (1981). "Western blotting": electrophoretic transfer of proteins from sodium dodecyl sulfate – polyacrylamide gels to unmodified nitrocellulose and radiographic detection with antibody and radioiodinated protein A. *Anal Biochem* **112**(2):195-203.
- Kurien, B.T. and Scofield, R.H. (2003). Protein blotting: a review. *J Imm Meth* **274**:1-15.
- Laemmli, U.K. (1970). Cleavage of structural proteins during the assembly of the head of bacteriophage T4. *Nature* **227**:680-5.
- Towbin, H., *et al.* (1979). Electrophoretic transfer of proteins from polyacrylamide gels to nitrocellulose sheets: procedure and some applications. *Proc Natl Acad Sci USA* **76**:4350-4.

This product ("Product") is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than the original purchaser of the Product ("Buyer").

No other warranties, express or implied, are granted, including without limitation, implied warranties of merchantability, fitness for any particular purpose, or non infringement. Buyer's exclusive remedy for non-conforming Products during the warranty period is limited to replacement of or refund for the non-conforming Product(s).

There is no obligation to replace Products as the result of (i) accident, disaster or event of force majeure, (ii) misuse, fault or negligence of or by Buyer, (iii) use of the Products in a manner for which they were not designed, or (iv) improper storage and handling of the Products.

Current product instructions are available at www.thermoscientific.com/pierce. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2014 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.