INSTRUCTIONS



PageRulerTM Unstained High Range Protein Ladder

Pub. No. MAN0011748 Rev. B.0 Pub Part No. 2162308.1

26637

Number Description

26637 PageRuler Unstained High Range Protein Ladder, 2 × 250μL

Storage Buffer: 62.5mM Tris•H₃PO₄ (pH 7.5 at 25°C), 1mM EDTA, 2% (w/v) SDS, 100mM DTT, 1mM NaN₃ 0.01% (w/v) bromophenol blue and 33% (v/v) glycerol.

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

Introduction

The Thermo Scientific PageRuler Unstained High Range Protein Ladder consists of a mixture of eight recombinant, purified proteins ranging from 60kDa to 250kDa. The ladder is visualized by SDS-PAGE using coomassie or silver stains or detected in Western blots with protein stains. For easy reference, the 150kDa protein band has a greater intensity than the other proteins in the ladder. 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.
- The large proteins (> 100kDa) in the ladder may require longer transfer times or higher transfer voltages for Western blotting.
- If additional bands appear in the protein ladder, add newly prepared dithiothreitol (DTT) solution to 100mM final concentration. DTT oxidation in the storage buffer can cause the appearance of additional bands.
- The amount of ladder can be reduced up to 10-fold for silver staining.

Procedure for Using the Protein Ladder in Polyacrylamide Gel Electrophoresis

- 1. Thaw the ladder at room temperature. Do not boil the protein ladder solution.
- 2. Mix the solution gently and thoroughly to ensure it 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)
 - Midi gel: 10μL per well (0.75-1.0mm thick) or 20μL per well (1.5mm thick)

Note: Dilute the ladder ~1/10 in reducing sample buffer for silver staining.

4. Return the unused protein ladder to -20°C for up to one year.



Related Products

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

26614 PageRuler Unstained Protein Ladder, 2 × 250µL 26616 PageRuler Prestained Protein Ladder, $2 \times 250 \mu L$ 26619 PageRuler Plus Prestained Protein Ladder, 2 × 250µL 26630 PageRuler Broad Range Unstained Protein Ladder, 2 × 250µL PageRuler Low Range Unstained Protein Ladder, 2 × 250µL 26632 26634 SpectraTM Multicolor Broad Range Protein Ladder, 2 × 250µL 26625 Spectra Multicolor High Range Protein Ladder, 2 × 250µL Spectra Multicolor Low Range Protein Ladder, 250µL 26628

LC5615 iBrightTM Prestained Protein Ladder

XP00060BOX NovexTM 6% Tris-Glycine Mini Gels, 10-well (see thermofisher.com/proteingels for a complete

listing)

EA0375BOX NuPAGETM 3-8% Tris-Acetate Protein Gels, 10-well (see thermofisher.com/proteingels for a

complete listing)

24615 Imperial™ Protein Stain, 1L LC6060 SimplyBlue™ SafeStain 24612 Pierce Silver Stain Kit

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 is manufactured under the license for Strep-tag® Technology covered by US patents 5,506,121 and 6,103,493 and foreign counterparts. Strep-tag is a trademark of the Institut fuer Bioanalytik GmbH.

For research use only. Not for use in diagnostic procedures.

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 thermofisher.com. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2017 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.