INSTRUCTIONS



PageRulerTM Prestained NIR Protein Ladder

Pub. No. MAN0011747 Rev. B.0 Pub Part No. 2162307.1

26635

Number

Description

26635 PageRuler Prestained NIR Protein Ladder, 2 × 250μL

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 NIR Protein Ladder is a mixture of 10 proteins ranging from 11kDa to 250kDa, which produces 10 well-defined blue bands by SDS-PAGE (see website for product images). The near-infrared (NIR) ladder can also be visualized by SDS-PAGE or Western blotting using a fluorescent detection instrument (e.g., TyphoonTM or LI-COR OdysseyTM). For easy reference, the 55kDa 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.
- Store the protein ladder for up to one year at -20°C.
- In low-percentage gels (< 10%), the low-molecular weight proteins in the ladder may migrate with the dye front.
- The large proteins (> 100kDa) 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 our website for migration patterns in different electrophoresis conditions.

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 (see Table 1).
- 4. Return the unused protein ladder to -20°C and store up to one year.
- 5. Fluorescently visualize the ladder using the 670nm red laser on a Typhoon Instrument or the 700nm channel on an Odyssey Instrument. (Optional)



Table 1. Volumes of Thermo Scientific PageRuler Prestained NIR Protein Ladder to load for different applications.

	Visual detection (0.75-1.0mm gel thickness)	Visual detection (1.5mm gel thickness)	Infrared* imaging detection	Western blotting detection
Mini-gel	5μL	10μL	1-5µL	1-3µL
Midi gel	10μL	$20\mu L$	2-10μL	2-6µL

^{*}For near-IR detection, optimize the ladder volume used.

Related Products

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

26614	PageRuler Unstained Protein Ladder, $2 \times 250 \mu L$
26616	PageRuler Prestained Protein Ladder, $2 \times 250 \mu L$
26619	PageRuler Plus Prestained Protein Ladder, $2 \times 250 \mu L$
26630	PageRuler Broad Range Unstained Protein Ladder, $2 \times 250 \mu L$
26632	PageRuler Low Range Unstained Protein Ladder, $2 \times 250 \mu L$
26634	Spectra TM Multicolor Broad Range Protein Ladder, $2 \times 250 \mu L$
26625	Spectra Multicolor High Range Protein Ladder, $2 \times 250 \mu L$
26628	Spectra Multicolor Low Range Protein Ladder, 250µL
LC5615	iBright™ Prestained Protein Ladder
XP04200BOX	Novex TM Tris-Glycine protein gels (see <u>thermofisher.com/proteingels</u> for a complete listing)
NW04120BOX	Bolt TM Bis-Tris Plus protein gels (see thermofisher.com/proteingels for a complete listing)

General References

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.

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.

Typhoon is a trademark of GE Healthcare Bio-sciences AB LLC.

Odyssey is a trademark of LI-COR Biosciences.

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.