

# PageRuler™ Low Range Unstained Protein Ladder

**26632**

Pub. No. MAN0011778

Rev. B.0

Pub Part No. 2162359.2

Number	Description
26632	PageRuler Low Range Unstained Protein Ladder, 2 × 250µL

**Storage Buffer:** 62.5mM Tris•H<sub>3</sub>PO<sub>4</sub> (pH 7.5 at 25°C), 1mM EDTA, 2% (w/v) SDS, 100mM DTT, 1mM NaN<sub>3</sub>, 0.01% (w/v) bromophenol blue and 33% (v/v) glycerol.

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

## Introduction

The Thermo Scientific PageRuler Low Range Unstained Protein Ladders are a mixture of seven recombinant proteins ranging from 5kDa to 100kDa and a synthetic peptide at 3.4kDa. This unstained ladder can be detected by coomassie or silver stains or in Western blots with protein stains. For easy reference, the 25kDa protein band is of greater intensity (see website for product images). The 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.
- In low-percentage gels (< 8%), low-molecular weight proteins (1-20kDa) may migrate with the dye front. Use high-percentage gels (>14%) to obtain well-resolved bands on low-molecular weight proteins
- Before coomassie or silver staining, fix proteins with 5% glutaraldehyde to preserve the low-molecular weight proteins in the ladder.
- For silver stain applications, dilute the protein ladder approximately 1/50 in reducing sample buffer.
- 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.

## Procedure for Use in Polyacrylamide Gel Electrophoresis

1. Thaw the protein ladder at room temperature. Do not boil ladder.
2. Mix the tube gently and thoroughly to ensure that the solution is homogeneous.
3. Load an appropriate volume of the protein 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)
4. Return the unused protein ladder to -20°C.

## Related Products

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

<b>26614</b>	<b>PageRuler Unstained Protein Ladder, 2 × 250µL</b>
<b>26616</b>	<b>PageRuler Prestained Protein Ladder, 2 × 250µL</b>
<b>26619</b>	<b>PageRuler Plus Prestained Protein Ladder, 2 × 250µL</b>
<b>26630</b>	<b>PageRuler Broad Range Unstained Protein Ladder, 2 × 250µL</b>
<b>26634</b>	<b>Spectra™ Multicolor Broad Range Protein Ladder, 2 × 250µL</b>
<b>26625</b>	<b>Spectra Multicolor High Range Protein Ladder, 2 × 250µL</b>
<b>26628</b>	<b>Spectra Multicolor Low Range Protein Ladder, 250µL</b>
<b>LC5615</b>	<b>iBright™ Prestained Protein Ladder</b>
<b>XP10200BOX</b>	<b>Novex™ 10-20% Tris-Glycine Mini Gels, 10-well</b> (see <a href="http://thermofisher.com/trisglycine">thermofisher.com/trisglycine</a> for a complete listing)
<b>NW04120BOX</b>	<b>Bolt™ 4-12% Bis-Tris Plus Gels, 10-well</b> (see <a href="http://thermofisher.com/bolt">thermofisher.com/bolt</a> for a complete listing)
<b>EC6625BOX</b>	<b>Novex 10-20% Tricine Protein Gels, 10-well</b> (see <a href="http://thermofisher.com/tricine">thermofisher.com/tricine</a> for a complete listing)
<b>24615</b>	<b>Imperial™ Protein Stain, 1L</b>
<b>LC6060</b>	<b>SimplyBlue™ SafeStain</b>

## 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.  
*Strep-tag*® is a trademark of IBA 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](http://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.