

Publication Part No. 10748010.pps Publication No. MAN0007901 Rev B.0

	<b>Package contents</b>	<b>Catalog Numbers</b> 10748-010	<b>Size:</b> 2 × 250 µL
---	-------------------------	-------------------------------------	----------------------------

	<b>Storage conditions</b>	<ul style="list-style-type: none"> <li>Store at -30°C to -10°C.</li> <li>Avoid repeated freezing and thawing.</li> </ul>
---	---------------------------	--

	<b>Required materials</b>	<ul style="list-style-type: none"> <li>Polyacrylamide gel(s)</li> <li>Electrophoresis apparatus for protein analysis</li> </ul>
---	---------------------------	---

	<b>Selection guide</b>	<p><a href="#">Pre-Stained Protein Standards</a> Go online to view related products.</p>
---	------------------------	--

- BenchMark™ Pre-Stained Protein Ladder allows you to accurately determine molecular weight and evaluate western transfer efficiency during electrophoresis.

- This standard consists of 10 pre-stained protein bands ranging in molecular weight from 6–180 kDa.

- Proteins are covalently coupled to blue dye, with one pink dyed band (4<sup>th</sup> from the top) to ensure correct identification of each protein band

- Storage buffer: 50 mM Tris-HCl (pH 6.8), 5 mM EDTA, 10 mM DTT, 1% (w/v) SDS, 10% (w/v) Glycerol.

- Use with Bolt™ Bis-Tris Plus, or Tris-Glycine gels. Click here to view our online [Protein Gel Selection Guide](#).

- DO NOT apply heat or add a reducing agent. BenchMark™ Pre-Stained Protein Ladder is supplied ready-to-use.



## Product description



## Important guidelines



## Online resources

Visit our [product page](#) for additional information and protocols. For support, visit [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).



For Research Use Only. Not for use in diagnostic procedures.

18 March 2014

## Protocol

- Thaw the standard at room temperature.
- Vortex gently to ensure the solution is homogeneous.
- Load the standard on the gel (see table for recommended volumes).

Gel Type	Application	
	Electrophoresis	Blotting
Mini gel	10 µL	5 µL
Midi gel	20 µL	10 µL

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.lifetechnologies.com/termsandconditions](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).

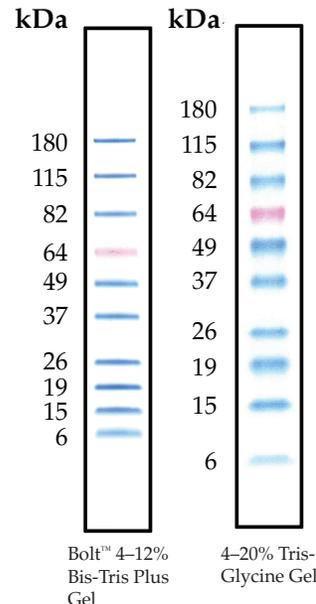
## Important licensing information

This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

## Disclaimer

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.



10 µL of BenchMark™ Pre-Stained Protein Ladder on gels with MES SDS Running Buffer



For support, visit [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).