

Publication Part No. 10747012.pps Publication No. MAN0007917 Rev B.0

 Package contents	Catalog Numbers 10747-012	Size: 2 × 250 µL
---	-------------------------------------	----------------------------

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

 Required materials	<ul style="list-style-type: none"> Polyacrylamide gel(s) Electrophoresis apparatus for protein analysis Appropriate sample buffer (when silver staining) 	
---	---	--

 Selection guide	Unstained Protein Standards Go online to view related products.	
--	--	--

- BenchMark™ Unstained Protein Ladder allows you to accurately determine molecular weight for polyacrylamide gel electrophoresis.
- This standard consists of 15 protein bands ranging in molecular weight from ~10–220 kDa.
- The 20 and 50 kDa protein bands are more prominent for easy orientation and identification.
- Storage buffer: 50 mM Tris-HCl (pH 6.8), 2 mM EDTA, 10 mM DTT, 2% (w/v) SDS, 10% (w/v) Glycerol, 0.01% (w/v) Bromophenol blue.

 Product description	<ul style="list-style-type: none"> The BenchMark™ Unstained Protein Ladder is supplied in gel loading buffer and can be applied directly to an SDS-polyacrylamide gel. Click here to view our online Protein Gel Selection Guide. 	
--	--	--

 Important guidelines	<ul style="list-style-type: none"> Visualize bands using Coomassie or silver staining techniques. DO NOT apply heat or add a reducing agent. BenchMark™ Unstained Protein Ladder is supplied ready-to-use. 	
---	--	--

 Online resources	Visit our product page for additional information and protocols. For support, visit 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.
- (*Silver staining only*) Dilute the standard 1:20 in 1X Sample Buffer.
- Load the standard on the gel (see table for recommended volumes).

Gel Type	Volume
Mini gel	5 µL
Midi gel	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. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

Important licensing information

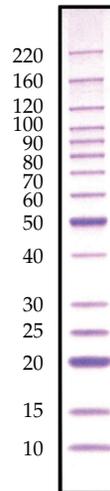
This products may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

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.

kDa



5 µL of BenchMark™ Protein Ladder on a NuPAGE® 4–12% Bis-Tris Gel with MES SDS Running Buffer, stained with Coomassie R-250.

life
technologies™

For support, visit www.lifetechnologies.com/support.