

### Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact [outlicensing@lifetech.com](mailto:outlicensing@lifetech.com) or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

©2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support) or email [techsupport@lifetech.com](mailto:techsupport@lifetech.com)

[www.lifetechnologies.com](http://www.lifetechnologies.com)



**novex**<sup>®</sup>  
by *life* technologies™

### BenchMark™ Pre-Stained Protein Ladder

**Cat. no.** 10748-010 **Size** 2 × 250 µL **Store at** -30°C to -10°C  
**Pub. Part no.** 10748010.pps **MAN0000876** **Rev. Date:** 14 October 2011

#### Description

The BenchMark™ Pre-Stained Protein Ladder is designed for use in SDS-PAGE using Tris-Glycine gels for estimating protein molecular weight, and evaluating western transfer efficiency. Covalently coupled dyes allow easy visualization of the ladder during, and after electrophoresis.

Features of the BenchMark™ Pre-Stained Protein Ladder:

- Consists of 10 pre-stained protein bands in the range of 6–180 kDa (proteins are covalently coupled to a blue or pink dye)
- Includes a protein coupled to a pink dye (fourth band from the top) to ensure proper identification of each protein band
- Supplied in a ready-to-use format

**Note:** Coupling of the chromophores to the proteins affects the apparent molecular weight of proteins in SDS-PAGE relative to unstained standards. Use pre-stained protein ladders to determine the approximate molecular weight of proteins. To accurately estimate the molecular weight of proteins, use unstained protein ladders.

#### Specifications

**Contents:** 2 × 250 µl of BenchMark™ Pre-Stained Protein Ladder  
**Storage Buffer:** 50 mM Tris-HCl, pH 6.8; 5 mM EDTA; 10 mM DTT; 1% (w/v) SDS; 10% (w/v) glycerol  
**Storage:** Store at -30°C to -10°C  
**Stability:** 1 year at -20°C

**Product Use** For research use only. Not for human or animal therapeutic or diagnostic use.

### Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact [outlicensing@lifetech.com](mailto:outlicensing@lifetech.com) or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

©2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support) or email [techsupport@lifetech.com](mailto:techsupport@lifetech.com)

[www.lifetechnologies.com](http://www.lifetechnologies.com)



**novex**<sup>®</sup>  
by *life* technologies™

### BenchMark™ Pre-Stained Protein Ladder

**Cat. no.** 10748-010 **Size** 2 × 250 µL **Store at** -30°C to -10°C  
**Pub. Part no.** 10748010.pps **MAN0000876** **Rev. Date:** 14 October 2011

#### Description

The BenchMark™ Pre-Stained Protein Ladder is designed for use in SDS-PAGE using Tris-Glycine gels for estimating protein molecular weight, and evaluating western transfer efficiency. Covalently coupled dyes allow easy visualization of the ladder during, and after electrophoresis.

Features of the BenchMark™ Pre-Stained Protein Ladder:

- Consists of 10 pre-stained protein bands in the range of 6–180 kDa (proteins are covalently coupled to a blue or pink dye)
- Includes a protein coupled to a pink dye (fourth band from the top) to ensure proper identification of each protein band
- Supplied in a ready-to-use format

**Note:** Coupling of the chromophores to the proteins affects the apparent molecular weight of proteins in SDS-PAGE relative to unstained standards. Use pre-stained protein ladders to determine the approximate molecular weight of proteins. To accurately estimate the molecular weight of proteins, use unstained protein ladders.

#### Specifications

**Contents:** 2 × 250 µl of BenchMark™ Pre-Stained Protein Ladder  
**Storage Buffer:** 50 mM Tris-HCl, pH 6.8; 5 mM EDTA; 10 mM DTT; 1% (w/v) SDS; 10% (w/v) glycerol  
**Storage:** Store at -30°C to -10°C  
**Stability:** 1 year at -20°C

**Product Use** For research use only. Not for human or animal therapeutic or diagnostic use.

### Directions

The BenchMark™ Pre-Stained Protein Ladder is supplied in a ready-to-use format.

**Do not heat** or add reducing agent to the ladder.

1. Thaw the ladder at room temperature.
2. Vortex gently to ensure the solution is homogenous.
3. Load the ladder into the well of an SDS gel using the loading volumes listed below to obtain the best results:

Gel Type	Electrophoresis	Blotting
Mini-Gel	10 µL	5 µL
Standard Gel	20 µL	10 µL

### Example

The average apparent molecular weights of the protein bands in the BenchMark™ Pre-Stained Protein Ladder are shown below. Ten microliters of the BenchMark™ Pre-Stained Protein Ladder was separated on a Novex® 4–20% Tris-Glycine gel.

	Band No.	Average Apparent Molecular Weight
1	1	~180 kDa
2	2	~115 kDa
3	3	~82 kDa
4	4	~64 kDa*
5	5	~49 kDa
6	6	~37 kDa
7	7	~26 kDa
8	8	~19 kDa
9	9	~15 kDa
10	10	~6 kDa

\*Orientation band (pink in color)

### Directions

The BenchMark™ Pre-Stained Protein Ladder is supplied in a ready-to-use format.

**Do not heat** or add reducing agent to the ladder.

1. Thaw the ladder at room temperature.
2. Vortex gently to ensure the solution is homogenous.
3. Load the ladder into the well of an SDS gel using the loading volumes listed below to obtain the best results:

Gel Type	Electrophoresis	Blotting
Mini-Gel	10 µL	5 µL
Standard Gel	20 µL	10 µL

### Example

The average apparent molecular weights of the protein bands in the BenchMark™ Pre-Stained Protein Ladder are shown below. Ten microliters of the BenchMark™ Pre-Stained Protein Ladder was separated on a Novex® 4–20% Tris-Glycine gel.

	Band No.	Average Apparent Molecular Weight
1	1	~180 kDa
2	2	~115 kDa
3	3	~82 kDa
4	4	~64 kDa*
5	5	~49 kDa
6	6	~37 kDa
7	7	~26 kDa
8	8	~19 kDa
9	9	~15 kDa
10	10	~6 kDa

\*Orientation band (pink in color)