thermo scientific

PRODUCT INFORMATION Thermo Scientific Spectra Multicolor High Range Protein Ladder

Pub. No. MAN0011775 Rev. Date 19 June 2019 (Rev. C.00)

Components	#26625	#26626	
Spectra Multicolor High Range Protein Ladder	2 x 250 µL	50 µL	

Store at -20 °C

www.thermofisher.com

www.thermoscientific.com/pierce

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

Introduction

The Thermo Scientific[™] Spectra[™] Multicolor High Range Protein Ladder is a prestained mixture of eight recombinant proteins ranging from 40 kDa to 300 kDa. Three different chromophores are bound to the proteins, producing a brightly colored ladder specifically designed for large protein analysis. The protein ladder is conveniently packaged and ready to use with no heating, diluting or additional reducing agent necessary.

Lot-to-lot variation of the apparent molecular weight of prestained proteins is \sim 5 %.

Storage Buffer: 62.5 mM Tris•H₃PO₄ (pH 7.5 at 25 °C), 1 mM EDTA, 2 % (w/v) SDS, 10 mM DTT, 1 mM NaN₃, 33 % (v/v) glycerol. Important Product Information

- Do not boil the protein ladder.
- The large proteins (> 100 kDa) in the protein ladder may require longer transfer times or higher transfer voltages for Western blotting.
- The mobility of proteins in the ladder 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 the table provided for migration patterns in different electrophoresis conditions.
- For precise MW determination use the Thermo Scientific[™] PageRuler[™] Unstained High Range Protein Ladder (#26637).

Gel type		Tris-Glycine						Tris-Acetate*		Bis-Tris*	
Gel concentration		4-12%	4%	6%	8%	10%	4-20%	3-8%	7%	4-12%	
Running buffer		Tris-Glycine						Tris-A	cetate	MOPS	
		Apparent Molecular Weights, kDa									
% lenght of gel	20	300 250 180 130 100 70 50 40	 300 250 180 130 100 	$ \begin{array}{c} 300 \\ 250 \\ 180 \\ - 130 \\ - 100 \\ - 70 \\ - 50 \\ - 40 \\ \end{array} $	 300 250 180 130 100 70 50 40 	300 250 180 130 - 100 - 70 - 50 - 40	 300 250 180 130 100 70 50 40 	 270 205 150 120 85 65 50 40 	 270 205 150 120 85 65 50 40 	 270 185 140 115 80 65 50 40 	

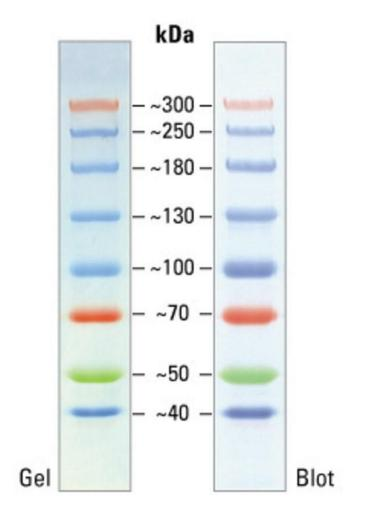
Migration Patterns of Spectra Multicolor High Range Protein Ladder

* migration patterns were determined using NuPAGE® precast gels.

Recommendations for Loading

- 1. Thaw the ladder at room temperature for a few minutes to dissolve precipitated solids. **Do not boil!**
- 2. Mix gently, but thoroughly, to ensure that the solution is homogeneous.
- 3. Load the following volumes of the ladder on an SDS-polyacrylamide gel:
 - 10 µL per well for mini gel,
 - $-\,20~\mu L$ per well for large gel.
 - Use the same volumes for Western blotting.
 - The loading volumes listed above are recommended for gels with a thickness of 0.75 mm-1.0 mm. The loading volume should be doubled for 1.5 mm thick gels.

Spectra Multicolor High Range Protein Ladder



4-12 % Tris-glycine SDS-PAGE

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.

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.

PRODUCT USE LIMITATION

This product is developed, designed and sold exclusively *for research purposes and in vitro use only.* The product was not tested for use in diagnostics or for drug development, nor is it suitable for administration to humans or animals. Please refer to <u>www.thermofisher.com</u> for Material Safety Data Sheet of the product.

 $\ensuremath{\mathbb{C}}$ 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.