

ProLite™ 5-245 kD Protein Ladder

Catalog number: 18015
Unit size: 100 uL

Component	Storage	Amount (Cat No. 18015)
ProLite™ 5-245 kD Protein Ladder	Freeze (< -15 °C), Minimize light exposure	100 µL

OVERVIEW

The ProLite™ 5-245 kD Protein Ladder is a high-quality three-color protein standard consisting of 13 prestained proteins. These proteins span a broad range of molecular weights, from 3.5 to 245 kDa. While most proteins in this ladder are covalently linked to a blue chromophore, it includes two distinctive reference bands – a green one at 25 kDa and a red one at 75 kDa that become evident when separated by SDS-PAGE. The ProLite™ 5-245 kD Protein Ladder is primarily used to track protein separation during SDS-polyacrylamide gel electrophoresis, confirm Western transfer efficiency across various membranes (like PVDF, nylon, or nitrocellulose), and to provide an estimation of protein sizes. It is conveniently supplied in a ready-to-use format premixed in a gel loading buffer. This protein ladder can be directly loaded onto gels without the need to heat, reduce, or introduce any reducing agents prior to its application.

AT A GLANCE

ProLite™ 5-245 kD Protein Ladder is stable for 12 months upon receiving when stored at <-15 °C with reduced light exposure.

SAMPLE EXPERIMENTAL PROTOCOL

Following usage protocol can be used as a guideline.

Sample protocol for usage in gels

1. Add 3 µL or 5 µL per loading for clear visualization during electrophoresis.
2. Add 1.5 µL to 2.5 µL per well for general western transferring on 15-well or 10-well mini-gel, respectively.
3. Apply more for thicker (>1.5 mm) or larger gel.

EXAMPLE DATA ANALYSIS AND FIGURES

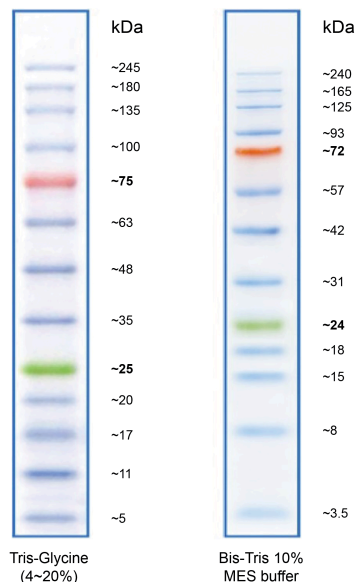


Figure 1. ProLite™ 5-245 kD Protein Ladder.

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.