Peacock™ Plus Prestained Protein Marker

A visible three-color protein marker for SDS-PAGE or western blotting, with 12 bands ranging from 8 to 245 kDa.

#Edition

Product attributes

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

PeacockTM Plus Prestained Protein Marker is a three-color protein ladder that allows you to visually monitor protein separation during SDS-PAGE or protein transfer to membranes for western blotting. We also offer PeacockTM Prestained Protein Marker, a three-color marker with 10 bands ranging from 10 kDa to 180 kDa

- 3-color visible protein ladder
- 12 total protein bands, ranging from 8 kDa to 245 kDa
- 10 blue bands, plus a red band at 75 kDa and green band at 25 kDa
- Ready-to-load: no heating required
- Store at -20°C

The PeacockTM Plus Prestained Marker contains a total of 12 visible bands. This includes 10 blue bands ranging from 8 kDa to 245 kDa, plus a red band at 75 kDa and green band at 25 kDa for easy band identification. PeacockTM Plus Prestained Protein Marker is ready to load with no heating or other preparation needed. Recommended loading is 3-5 uL per well for mini-gels.

Also visit our <u>Protein Detection</u>, <u>Quantitation</u>, <u>& Analysis</u> technology page to learn about our safe and sensitive protein gel stains and western blotting normalization reagents.

This datasheet was generated on November 3, 2025 at 04:06:22 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/peacock-plus-prestained-protein-marker/