



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## Peacock™ Prestained Protein Marker

A visible three-color protein marker for SDS-PAGE or western blotting, with 10 bands ranging from 10 to 180 kDa.



### Product attributes

## Product Description

Peacock™ 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 [Peacock™ Plus Prestained Protein Marker](#), a three color marker with 12 bands ranging from 8 kDa to 245 kDa.

- 3-color visible protein ladder
- 10 total protein bands, ranging from 10 kDa to 180 kDa
- 8 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 Peacock™ Prestained Marker contains a total of 10 visible bands. This includes 8 blue bands ranging from 10 kDa to 180 kDa, plus a red band at 75 kDa and green band at 25 kDa for easy band identification. Peacock™ 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 [Protein Detection, Quantitation, & Analysis](#) technology page to learn about our safe and sensitive protein gel stains and western blotting normalization reagents.

This datasheet was generated on January 8, 2026 at 03:18:25 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/peacock-prestained-protein-marker/>