

# Product Information

## Peacock™ Plus Prestained Protein Marker

**Catalog Number:** 21531-50uL, 21531-500uL

**Unit Size:** 50 uL or 500 uL

### Storage and Handling

Store at -20°C. Product is stable for at least one year from date of receipt when stored as recommended. Alternatively, the marker can be stored at room temperature for up to 2 weeks, or at 4°C for up to three months.

### Product Description

Peacock™ Plus Prestained Protein Marker is a three-color protein ladder. It allows visual monitoring of protein separation during SDS-PAGE and protein transfer to membranes for western blotting, as well as estimation of protein molecular weight. It contains 10 visible blue bands ranging from 8 to 245 kDa, plus red and green bands at 75 kDa and 25 kDa for easy band identification.

The marker is provided in 10 mM Tris-phosphate buffer (pH 7.5 at 25°C), 2% SDS, 0.2 mM DTT, 3.6 M Urea, and 15% glycerol. Warm to room temperature and mix before use. Do not heat, dilute, or add reducing agent before loading. The recommended loading volume is 3-5 uL per well for mini-gels.

We also offer Peacock™ Prestained Protein Marker, with 10 bands ranging from 10-180 kDa. Also see our rapid, easy-to-use, and non-toxic One-Step Blue™ and One-Step Lumitein™ protein gel stains (see related products).

### Related Products

Catalog No.	Product
21530	Peacock™ Prestained Protein Marker
21003	One-Step Blue™ Protein Gel Stain
21004	One-Step Lumitein™ Protein Gel Stain
21005	One-Step Lumitein™ UV Protein Gel Stain
22001	Ponceau S Solution
30071	AccuOrange™ Protein Quantitation Kit
22010	10X Fish Gelatin Blocking Agent
22012	Non-fat dry milk
22011	Fish gelatin powder
22014	BSA, IgG- and protease-free, 30% solution
22002	TWEEN® 20
41003	GelRed™ Nucleic Acid Gel Stain
41005	GelGreen™ Nucleic Acid Gel Stain
41008	PAGE GelRed™ Nucleic Acid Gel Stain

Please visit our website at [www.biotium.com](http://www.biotium.com) for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, fluorescent CF™ dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.

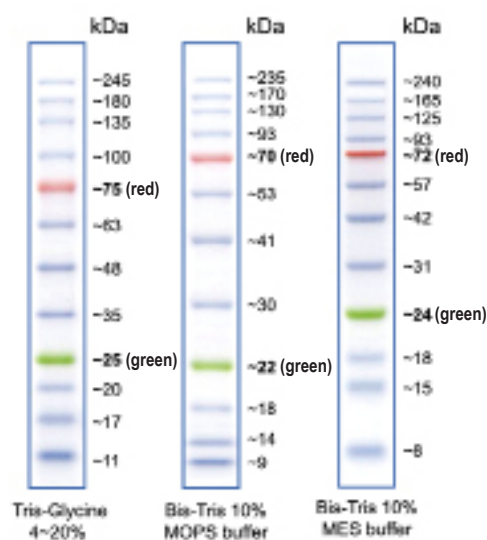


Figure 1. Apparent molecular weights of bands in Peacock™ Plus Prestained Protein Marker on various SDS-PAGE gels.