Revised: September 21, 2022

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 orange and green bands at 75 kDa and 25 kDa, respectively, for easy band identification (Figure 1).

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.

Peacock™ Plus Prestained Protein Marker can be used with near-infrared (NIR) fluorescent western blots. NIR fluorescence will be readily detectable for the blue and green marker bands in the 700 nm detection channel, while the orange marker band will be faint or undetectable (Figure 2). Marker bands may be faintly detectable in the 800 nm detection channel at high gain/intensity settings. The amount of marker loaded on the gel may need to be adjusted to match the brightness of the antibody signal. If antibody signal is very bright for your samples, we recommend loading 3 uL of ladder to reach the appropriate brightness on NIR scanners. If your antibody signal is dimmer, we recommend loading 1.5 uL of ladder, so that the ladder signal does not overwhelm sample signal.

We also offer Peacock™ Prestained Protein Marker, with 10 bands ranging from 10-180 kDa. In addition, see our rapid, easy-to-use, and non-toxic One-Step protein gel stains (see Related Products).

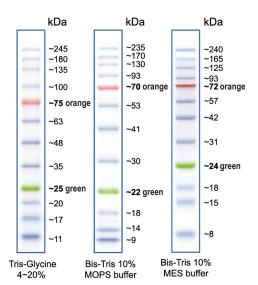


Figure 1. Apparent molecular weights of bands in Peacock™ Plus Prestained Protein Marker on various SDS-PAGE gels.

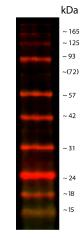


Figure 2. Peacock™ Plus Prestained Protein Marker, 3 uL loaded, after transfer to a western blot membrane. Imaged on a LI-COR Odyssey® infrared imaging system in the 700 nm channel. The blue and green marker bands are detectable, but the orange 72 kDa band is faint or undetectable by NIR fluorescence.

Related Products

Cat. No. Product	
	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
23013	TrueBlack® WB Blocking Buffer Kit
22010	10X Fish Gelatin Blocking Agent
22012	Dry Milk Powder
22011	Fish Gelatin Powder
22014	Bovine Serum Albumin 30% Solution
22002	TWEEN® 20
20192	CF®680R Goat Anti-Mouse IgG (H+L), Highly Cross-Adsorbed
20193	CF®680R Goat Anti-Rabbit IgG (H+L), Highly Cross-Adsorbed
20078	CF®770 Goat Anti-Rabbit IgG (H+L), Highly Cross-Adsorbed
20077	CF®770 Goat Anti-Mouse IgG (H+L), Highly Cross-Adsorbed
29072	CF®680R Streptavidin Conjugate
30071	AccuOrange™ Protein Quantitation Kit
33025	VersaBlot™ Total Protein Normalization Kits
40136	4X Protein Loading Buffer with Orange Tracking Dye

Please visit our website at 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.

TWEEN® is a registered trademark of Croda International LLC; Odyssey is a registered trademark of LI-COR, Inc.

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