

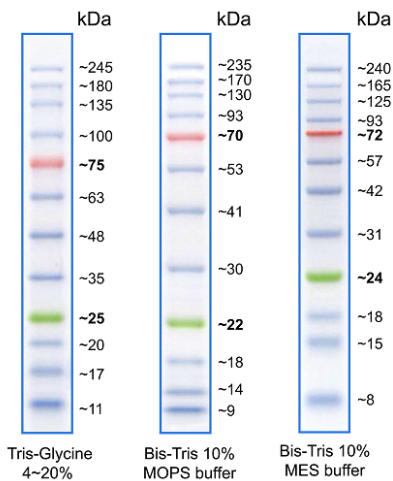


BlueEye Prestained Protein Marker

10-245 kDa

Ready-to-use prestained protein ladder

Cat. No.	Amount
PS-104	2 x 250 µl



Applications:

- Visualization of protein migration during SDS-polyacrylamide gel electrophoresis
- Verification of protein transfer onto membranes after Western blotting
- Approximate sizing of proteins on SDS-polyacrylamide gels and Western blots

Description:

The BlueEye Prestained Protein Marker is a three-color protein standard with 12 pre-stained purified proteins. Proteins are covalently coupled with a blue chromophore except for two reference bands, one green at 25 kDa and one red one at 75 kDa respectively, when separated on SDS-PAGE (Tris-glycine buffer).

Recommended Load:

3-5 µl Protein Molecular Weight Marker per lane

Migration Patterns

For *in vitro* use only!

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term storage (up to 1 month) at 4 °C possible.

Shelf Life: 12 months

Purity: > 95 % (SDS-PAGE)

Form: liquid (Supplied in 20 mM Tris-phosphate pH 7.5 , 2 % SDS, 0.2 DTT, 3.6 M urea and 15 % glycerol)