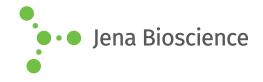
## **DATA SHEET**

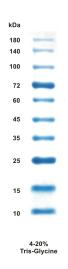




## ■ BlueClassic Prestained Protein Marker (10-180kDa)

10 - 180 kDa prestained protein ladder Ready-to-use protein ladder

Cat. No.	Amount
PS-107	2 x 250 μl



Migration Pattern

For general laboratory use.

**Shipping:** shipped on gel packs

Storage Conditions: store at -20  $^{\circ}\text{C}$ 

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  $\,$ 

mM DTT, 3.6 M urea and 15 % glycerol)

## **Description:**

The BlueClassic Prestained Protein Marker is a blue-colored protein standard with 10 pre-stained and purified proteins supplied in gel loading buffer for direct application in SDS-polyacrylamide gel.

## **Recommended Load:**

 $3-5~\mu l$  Protein Molecular Weight Marker per lane