

BSA (10 mg/ml)

Catalog No: 37497

Format: 0.1 ml

Quality Control: Active Motif's BSA (part number 104892), also known as bovine serum albumin, is a globular protein used to reduce non-specific binding events. BSA can also be used to stabilize some proteins.

BSA is supplied at a concentration of 10 mg/ml in 10 mM Tris-HCl, 50 mM KCl, 0.05 mM EDTA and 25% glycerol.

Storage and Guarantee: Store at -20°C. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.