



Product Information Sheet

Product Name: BSA-Biotin Conjugate

Catalog Number: AS-72264

Amount: 10 mg

Degree of Substitution: See label on the vial.
(DOS)

Formulation: 10 mM phosphate, 150 mM sodium chloride, pH 7.2.

Storage: This product is supplied as lyophilized powder and is stable for one year at -20°C. Solutions can be made by dissolving the conjugate powder in the sterile water. Store the solutions at 4°C with the addition of sodium azide at a final concentration of 0.01%. For long-term storage, divide the solution into aliquots and store at -20°C or add an equal volume of glycerol (ACS grade or higher) and store at -20°C without aliquoting. Avoid multiple thaw-freeze cycles.

Description: Bovine serum albumin (BSA) is an albumin protein with molecular weight of 66 kDa. Biotinylated BSA is widely used for numerous applications such as, binding assays¹⁻³, protein printing⁴, carrier protein test⁵ and as a permeability probe⁶.

References:

1. Hetland O., Scand J Clin Lab Invest. **55**(1995):701-713.
2. Emans N., Biophys J. **69**(1995):716-728.
3. Wu FB., Clin Chim Acta. **308**(2001):117-126.
4. Wilson WC Jr., Anat Rec A Discov Mol Cell Evol Biol. **272**(2003):491-496.
5. Lövgren K., J Immunol Methods. **98**(1987):137-143.
6. Tong H., Inflammation. **37**(2014): 27-34

For in vitro research use only.