

BSA

Albumin from bovine serum $\geq 98\%$

Cat.-No.	Amount
CPP-A02	1 g

For *in vitro* use only
Quality guaranteed for 12 months
Store at 4°C

Product Information

CAS Number: 9048-46-8

Solubility

soluble in water

Molecular Weight

66.4 kDa

BSA is a single polypeptide chain which consist of 583 amino acid residues and no carbohydrates. It contains 17 intrachain disulfide bridges and 1 sulfhydryl group at pH 5-7.

Description

Albumins are frequently used as stabilizers for other solubilized proteins (e.g., labile enzymes). BSA can act as co-substrate.

It reduces the proteolytic degradation of CPPs. We recommend concentrations of 0,5 to 1% BSA in serum free medium. BSA can stimulate the cargo internalisation, especially in tumor cells.

For more information see **general manual (4.2.2)**.

Selected References

The Plasma Proteins: Structure, Function and Genetic Control, 2nd, Vol. 1, p. 141, 147, Academic Press, New York (1975).
Reed, R.G. *et al.* (1980) Sequence of residues 400-403 of bovine serum albumin. *Biochem. J.*, **191**:867
Hirayama, K. *et al.* (1990) Rapid confirmation of the primary structure of bovine serum albumin by ESIMS and FRIT-FAB LC/MS. *BBRC*, **173**:639