

Atto488-BSA

Bovine serum albumin fluorescent-labeled with Atto488

Cat.-No.	Amount
CPP-A08S	100 µg
CPP-A08L	500 µg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C protected from light

Product Information

Bovine serum albumin (BSA) labeled by coupling with a sixfold molar excess of Atto488 N-hydroxy-succinimide ester.

Molecular Weight_{approx.}

68 kDa

λ_{exc}

501 nm

λ_{em}

523 nm

Solubility

soluble in water (1 µg/1 µl)

Description

Dissolve 100 µg in 100 µl sterile and oxygen free water, keep aliquots at -20°C. The required amount of Atto488-BSA depends on the aim of your experiments and on the sensitivity of your measurements. We recommend starting with 1 µg.

For complex formation, transduction, improving transduction and detection read **general manual**.