

**Atto488-BSA**

Bovine serum albumin fluorescent-labeled with Atto488

Cat. No.	Amount
CPP-A08L	500 µg

For in vitro use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** protect from light**Shelf Life:** 12 months after date of delivery**Molecular Weight:** 68 kDa**Form:** lyophilised orange powder**Solubility:** soluble in water (1 mg/ml)**Spectroscopic Properties:** λ_{exc} 501 nm; λ_{em} 523 nm;
 ϵ 90.0 L mmol⁻¹ cm⁻¹**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**.

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

Exclusively distributed in Japan by Greiner BioOne