



Product Information Sheet

Product Name: BSA-HiLyte™ Fluor 488 Conjugate

Catalog Number: AS-72265

Amount: 5 mg

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

Spectral properties: Green fluorescence. Excitation/Emission wavelength= 499 nm/523 nm

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 longer storage, make the aliquots and store at -20°C. Avoid multiple thaw-freeze cycle and protect from light.

Description: Bovine serum albumin (BSA) fluorescent conjugates are commonly used in various applications such as microvascular imaging¹, electroporation quantification², receptor-mediated endocytosis/exocytosis³ and drug delivery studies⁴.

The fluorescence of HiLyte™ Fluor 488-BSA conjugate can be observed at the excitation/emission wavelength of 499 nm/523 nm.

References:

1. Lugo-Hernandez E., J Cereb Blood Flow Metab. **37**(2017):3355-67.
2. Muñoz CF., Algal Research. **33**(2018):448-55.
3. Weaver DJ Jr., Biol Cell. **90**(1998):169-81.
4. Wang Q., J Pharm Sci. **100**(2011):4804-14.

For in vitro research use only.