



Product Information Sheet

- Product Name:** BSA-HiLyte™ Fluor 647 Conjugate
- Catalog Number:** AS-72267
- Amount:** 5 mg
- Degree of Substitution: (DOS)** See label on the vial
- Spectral properties:** Far-red fluorescence. Excitation/Emission wavelength= 649 nm/672 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 647-BSA conjugate can be observed at the excitation/emission wavelength of 649 nm/672 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. Campana PT., Bull Chem Soc Jpn. **92**(2019):1237-44.

For *in vitro* research use only.