



## Product Information Sheet

---

- Product Name:** BSA-HiLyte™ Fluor 555 Conjugate
- Catalog Number:** AS-72266
- Amount:** 5 mg
- Degree of Substitution: (DOS)** See label on the vial
- Spectral properties:** Orange fluorescence. Excitation/Emission wavelength= 550 nm/566 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 imaging<sup>1</sup>, electroporation quantification<sup>2</sup>, receptor-mediated endocytosis/exocytosis<sup>3</sup> and drug delivery studies<sup>4</sup>.
- The fluorescence of HiLyte™ Fluor 555-BSA conjugate can be observed at the excitation/emission wavelength of 551 nm/567 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.**