



## Product Information Sheet

---

<b>Product Name:</b>	BSA-HiLyte™ Fluor 750 Conjugate
<b>Catalog Number:</b>	AS-72268
<b>Amount:</b>	5 mg
<b>Degree of Substitution: (DOS)</b>	See label on the vial
<b>Spectral properties:</b>	Near-infrared fluorescence. Excitation/Emission wavelength= 754 nm/778 nm
<b>Formulation:</b>	10 mM phosphate, 150 mM sodium chloride, pH 7.2.
<b>Storage:</b>	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.
<b>Description:</b>	<p>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>.</p> <p>The fluorescence signal of HiLyte™ Fluor 750-BSA conjugate can be observed at the excitation/emission wavelength of 754 nm/778 nm.</p>

### 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.**