



## Product Information Sheet

---

<b>Product Name:</b>	OVA-HiLyte™ Fluor 488 Conjugate
<b>Catalog Number:</b>	AS-72270
<b>Amount:</b>	1 mg
<b>Degree of Substitution: (DOS)</b>	See label on the vial
<b>Spectral properties:</b>	Green Fluorescence. Excitation/Emission wavelength= 499 nm/523 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>Ovalbumin (OVA) is the major protein found in the egg white. The fluorescent dye conjugated OVA have been used in antigen uptake assay <sup>1,2</sup>, lymphatic circulation tracking <sup>3</sup>and drug delivery studies<sup>4</sup>.</p> <p>The fluorescence of HiLyte™ Fluor 488-OVA conjugate can be observed at the excitation/emission wavelength of 499 nm/523 nm.</p>
<b>References:</b>	<ol style="list-style-type: none"><li>1. Garrett WS., Cell. <b>10</b>(2000):325-34.</li><li>2. Granucci F., Microbes Infect. <b>1</b>(1999):1079-84.</li><li>3. Lin Y., J Immunol. <b>201</b>(2018):296-305.</li><li>4. Akagi T., Biomaterials. <b>28</b>(2007):3427-36.</li></ol>

**For *in vitro* research use only.**