



Product Information Sheet

- Product Name:** OVA-HiLyte™ Fluor 647 Conjugate
- Catalog Number:** AS-72272
- Amount:** 1 mg
- Degree of Substitution: (DOS)** See label on the vial
- Spectral properties:** Far-red fluorescence. Excitation/Emission wavelength= 650 nm/675 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:** Ovalbumin (OVA) is the major protein found in the egg white. The fluorescent dye conjugated OVA have been used in antigen uptake assay^{1,2}, lymphatic circulation tracking³ and drug delivery studies⁴.

The fluorescence of HiLyte™ Fluor 647-OVA conjugate can be observed at the excitation/emission wavelength of 650 nm/675 nm.
- References:**
1. Garrett WS., Cell. **10**(2000):325-34.
 2. Granucci F., Microbes Infect. **1**(1999):1079-84.
 3. Lin Y., J Immunol. **201**(2018):296-305.
 4. Akagi T., Biomaterials. **28**(2007):3427-36.

For *in vitro* research use only.