



Product Information Sheet

Product Name: OVA-Biotin Conjugate

Catalog Number: AS-72269

Amount: 10 mg

Degree of Substitution: See label on the vial.
(DOS)

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 long-term storage, divide the solution into aliquots and store at -20°C or add an equal volume of glycerol (ACS grade or higher) and store at -20°C without aliquoting. Avoid multiple thaw-freeze cycles.

Description: Ovalbumin (OVA) is the major protein found in the egg white. The biotinylated OVA have been used in binding assays¹⁻³, immune complex recycling assays⁴, and as a protein analysis standard⁵.

References:

1. Arribillaga L., Biomed Res Int. 2013:864720.
2. Heron BT., Int J Parasitol. **41**(2011):137-144.
3. Cao Y., J Immunol Methods. **220**(1998):85-91.
4. Bergtold A., Immunity. **23**(2005):503-514.
5. O'Hara JM., J Immunol Methods. **395**(2013):71-78.

For *in vitro* research use only.