## **cOvalbumin: Cationized with Ethylenediamine**

Catalog Number:	5605
Size:	10 mg
Molecular Weight:	44 kD
Solubility:	Soluble in water, giving a clear or yellowish solution.
Description:	cOvalbumin is prepared by treating the native ovalbumin with ethylenediamines that replace most of negatively-charged carboxyl groups with positively-charged primary amines, resulting a highly positively-charged cOvalbumin. The cationization significantly increases the immunogenicity compared to native ovalbumin. In addition, the increased number of primary amines provides more conjugation sites available for hapten molecules with general conjugation methods.
Storage/Stability:	Store at -20°C/1 year
Format:	Lyophilized in PBS, pH 7.2.
Immunogen:	Use as a carrier protein for immunization
Purification:	cOvalbumin is purified by a fractionation method, and is supplied with the purity over 97% by SDS.
Applications:	cOvalbumin itself acts an excellent immnogen with a greater immunogenicity compared to the native ovalbumin. With increased number of free amines, more antigen molecules can be coupled to cOvalbumin. When a stronger immunogenicity and a high concentration of hapten are needed, cOvalbumin is a good choice for the immunogen preparation of small hapten molecules.
References:	Sheng-Liang Deng, Ping Li, Hong-Bin Liu, Shu-Ming Yang (2014) Preparation and characterization of ultrasensitive and specific polyclonal antiserum against ciprofloxacin based on cationized bovine serum albumin. <b>Chemical Papers</b> 68 (11) 1505–1513.