

Ovalbumin

Chicken egg albumin

Catalog # vac-stova

<http://www.invivogen.com/ovalbumin>

For research use only.

Version # 17E02-MM

PRODUCT INFORMATION

Content:

- 1 g ovalbumin

Storage and stability

- Ovalbumin is shipped at room temperature and should be stored at 4°C.
- Upon resuspension, prepare aliquots of ovalbumin and store at -20°C. Resuspended product is stable for 6 months when properly stored. Avoid repeated freeze-thaw cycles.

Quality control

- ≥ 95% purity (agarose gel electrophoresis)
- Functionality of ovalbumin was assessed by enzyme-linked immunosorbent assay (ELISA).
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

DESCRIPTION

Ovalbumin (also known as albumin or OVA) is a key reference protein for immunization and biochemical studies. Ovalbumin is the major protein constituent of chicken egg whites. Ovalbumin is a phosphorylated-glycoprotein.

RELATED PRODUCTS

Product	Description	Catalog Code
Alum and Emulsions		
AddaVax™	Squalene-Oil-in-water	vac-adx-10
Alhydrogel 2%	Aluminium hydroxide gel	vac-alu-250
IFA	Incomplete Freund's adjuvant	vac-ifa-10
PRR Ligands		
3'3'-cGAMP VacciGrade™	STING agonist	vac-nacga
Flagellin FlIC VacciGrade™	TLR5 agonist	vac-fla
Gardiquimod VacciGrade™	TLR7 agonist	vac-gdq
MPLA VacciGrade™	TLR4 agonist	vac-mpl
N-glycyl-MDP VacciGrade™	NOD2 agonist	vac-gmdp
ODN 1585 VacciGrade™	murine TLR9 agonist	vac-1585-1
Pam3CSK4 VacciGrade™	TLR2 agonist	vac-pms
Poly(I:C) VacciGrade™	TLR3 agonist	vac-pic
R848 VacciGrade™	TLR7/8 agonist	vac-r848
OVA Antigens		
Ova 257-264	For detection; ELISPOT	vac-sin
Ova 323-339	For detection; ELISPOT	vac-isq

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873
InvivoGen USA (International): +1 (858) 457-5873
InvivoGen Europe: +33 (0) 5-62-71-69-39
InvivoGen Hong Kong: +852 3-622-34-80
E-mail: info@invivogen.com

METHODS

Preparation of stock solution (10 mg/ml)

- Weigh 10 mg of ovalbumin.
- Add 1 ml of **room-temperature** water to 10 mg of ovalbumin to obtain a solution at 10 mg/ml.
- Mix the solution by pipetting up and down. The solution may appear slightly hazy. Further dilutions can be made with water or phosphate buffered saline (PBS).

Note: Allow water or PBS to reach room temperature before use.

CHEMICAL PROPERTIES

CAS number: 9006-59-1

Molecular weight: ~ 45 kDa

Solubility: 40 mg/ml in water

Appearance: White to faint yellow

APPLICATIONS

Ovalbumin can be used for ELISA at a working concentration of 10 - 50 µg/ml in PBS (pH 7.4).

Other applications have not been tested.