

OVA 257-264

Chicken egg albumin peptide

Catalog code: vac-sin

<https://www.invivogen.com/ova-peptide>

For research use only

Version 22F27-MM

PRODUCT INFORMATION

Contents

- 1 mg OVA 257-264 provided lyophilized

Storage and stability

- OVA 257-264 is shipped at room temperature. Upon receipt, store at -20°C.

Note: OVA 257-264 is hygroscopic and light sensitive. Keep tightly closed and protect from light.

- Upon resuspension, prepare aliquots of OVA 257-264 and store at -20°C. Resuspended product is stable for 6 months when properly stored. Avoid repeated freeze-thaw cycles.

Quality Control

- Purity: ≥ 95% (UHPLC)
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

DESCRIPTION

OVA 257-264 is a class I (Kb)-restricted peptide epitope of ovalbumin (OVA), presented by the class I major histocompatibility complex (MHC) molecule, H-2Kb¹. It has been demonstrated that OVA 257-264 peptides can be used to detect a strong CD8+ cytolytic T cell response².

Sequence:

SIINFEKL (H-Ser - Ile - Ile - Asn - Phe - Glu - Lys - Leu - OH)

1. Rötzschke O. *et al.*, 1991. Exact prediction of a natural T cell epitope. *Eur J Immunol.* 21(11):2891-4. 2. Lipford G.B. *et al.*, 1993. Primary in vivo responses to ovalbumin. Probing the predictive value of the Kb binding motif. *J Immunol.* 150(4):1212-1222.

CHEMICAL PROPERTIES

CAS number: 138831-86-4

Synonym: SIINFEKL

Chemical formula: C₄₅H₇₄N₁₀O₁₃

Molecular weight: 963.2 g/mol

Solubility: 1 mg/ml in water

METHODS

Preparation of stock solution (1 mg/ml)

1. Add 1 ml water to 1 mg of OVA 257-264.
2. Mix the solution by pipetting up and down.

APPLICATIONS

OVA 257-264 can be used for Enzyme-Linked ImmunoSpot (ELISPOT) assay at a working concentration of 10 µg/ml.

Other applications have not been tested.

RELATED PRODUCTS

Product	Description	Cat.Code
EndoFit™ Ovalbumin	Ova for <i>in vivo</i> use	vac-pova
Ovabumin	Ova for detection use	vac-stova
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 Asia: +852 3622-3480

E-mail: info@invivogen.com