

## **Product Data Sheet**

Product Name: OVA (323-339), amide

Catalog Number: AS-62571-1 (1 mg) Lot Number: See label on vial

AS-62571-5 (5 mg)

Sequence: H-Ile-Ser-Gln-Ala-Val-His-Ala-Ala-His-Ala-Glu-Ile-Asn-Glu-Ala-Gly-Arg-

NH2 (3-letter code)

ISQAVHAAHAEINEAGR-NH2 (1-letter code)

Molecular Weight: 1774.0

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: OVA (323-339) peptide is freely soluble in H<sub>2</sub>O.

Storage: OVA (323-339) peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at –20°C or lower. Reconstituted peptide can be aliquoted and stored at –20°C or lower.

Description: This peptide is amino acids 323 to 339 amidated fragment of ovalbumin (OVA), the H-2b-restricted OVA class II epitope. This peptide is recognized by many T cells because it contains multiple T cell epitopes. This OVA fragment contains a nested set of CD4+ T cell epitopes. OVA 323 to 339 can be presented by I-Ad in at least three binding registers. The residues 327 to 333 are critical for peptide binding to I-Ad. Ref: Robertson, J. et al. *J. Immunol.* **164**, 4706 (2000), Oran, A. and Robinson, H. *J. Immunol.* **171**, 1999 (2003), Nag, B. et al. *J. Immunol.* **148**, 2040 (1992), Hongming, Z. et al. *J. Bacteriol.* **187**, 175 (2005).

## Related Products:

Name	Cat #	Size
OVA (323-339)	AS-27024	1 mg
(ISQAVHAAHAEINEAGR)	AS-27025	5 mg
OVA (257-264)	AS-60193-1	1 mg
(SIINFEKL)	AS-60193-5	5 mg

For Research Use Only