

PRODUCT INFORMATION



Ovalbumin (257-264) (chicken) (trifluoroacetate salt)

Item No. 36882

CAS Registry No.: 1262751-08-5

Formal Name: L-seryl-L-isoleucyl-L-isoleucyl-L-asparaginyl-L-phenylalanyl-L-alpha-glutamyl-L-lysyl-L-leucine, monotrifluoroacetate salt

Synonyms: Ser-Ile-Ile-Asn-Phe-Glu-Lys-Leu, SIINFEKL, OVA (257-264)

MF: $C_{45}H_{74}N_{10}O_{13} \cdot CF_3COOH$

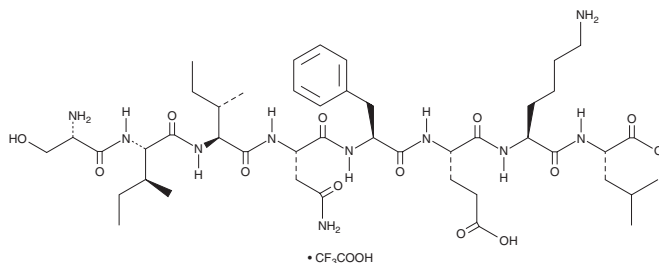
FW: 1,007.2

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Ovalbumin (257-264) (chicken) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the ovalbumin (257-264) (chicken) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Ovalbumin (257-264) (chicken) (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ovalbumin (257-264) (chicken) (trifluoroacetate salt) in these solvents is approximately 2 mg/ml.

Description

Ovalbumin (257-264) is a synthetic peptide fragment corresponding to the mouse MHC class I (H-2K^b) CD8⁺ T cell-restricted epitope of chicken ovalbumin.^{1,2} It is presented by H-2K^b dominant T cells and has been used to induce CD8⁺ T cell responses *in vitro* and *in vivo*.¹⁻³

References

1. Wick, M.J. and Pfeifer, J.D. Major histocompatibility complex class I presentation of ovalbumin peptide 257-264 from exogenous sources: Protein context influences the degree of TAP-independent presentation. *Eur. J. Immunol.* **26**(11), 2790-2799 (1996).
2. Chaoul, N. and Dadglio, G. Antigen-specific *in vivo* killing assay. *Methods Mol. Biol.* **2325**, 55-64 (2021).
3. Berguer, P.M., Alzogaray, V.A., Rossi, A.H., et al. A polymeric protein induces specific cytotoxicity in a TLR4 dependent manner in the absence of adjuvants. *PLoS One* **7**(9), e45705 (2012).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/21/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM