

Product Data Sheet

Product Name: Influenza A NP(366-374) Strain A/PR/8/35

Catalog Number: AS-60624 (0.5 mg) Lot Number: See label on vial

Sequence: H-Ala-Ser-Asn-Glu-Asn-Met-Glu-Thr-Met-OH (3-letter code)

ASNENMETM (1-letter code)

Molecular Weight: 1027.1

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Influenza A NP peptide is freely soluble in water.

Storage: Influenza A NP 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 an H2-Db-restricted epitope from the Influenza A/PR/8/34 nucleoprotein. Ref: Mo, X. et al. *J. Immunol.* **163**, 5851 (1999); Haanen, J. et al. *J. Exp. Med.* **90**, 1319 (1999); Bennett N. et al. *PLoS. Biol.* **3**, e120 (2005).

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

RSV M2₈₂₋₉₀ (SYIGSINNI) and RSV M₁₈₇₋₁₉₅ (NAITNAKII) were derived from the M2 and M proteins, respectively, of the RSV A2 strain. An H2-K^d-binding influenza virus A/Puerto Rico/8/34 nucleoprotein peptide comprised of residues 147 to 155 (NP₁₄₇₋₁₅₅; TYQRTRALV) and an H2-D^b-binding influenza A/Puerto Rico/8/34 peptide NP₃₆₆₋₃₇₄ (ASNENMETM) were used as negative controls. All peptides were synthesized by Anaspec, Inc. (San Jose, CA), and confirmed to be >95% pure by analytical high-performance liquid chromatography at the National Institute of Allergy and Infections Diseases peptide core facility (Bethesda, MD). For negative controls, cells from uninfected mice were stained, and phycoerythrin-labeled tetrameric complexes of H-2D^b with the influenza virus NP₃₆₆₋₃₇₄ peptide, H-2K^d with the influenza virus NP₁₄₇₋₁₅₅ peptide, or H-2K^b with an ovalbumin epitope comprised of residues 257 to 264 (Ova₂₅₇₋₂₆₄; SIINFEKL) were used-Ruckwardt, T. et al. J. *Virol.* 10.1128/JVI.00036-09 (2009).

Published Citations:

Ruckwardt, T. et al. J. Virol. 10.1128/JVI.00036-09 (2009).

Related Products:

Name	Cat #	Size
Influenza A NP (366-374) Strain A/NT/60/68	AS-60623	1 mg
ASNENMDAM		

For Research Use Only