



Product Data Sheet

Product Name:	HA peptide	
Catalog Number:	AS-21156 (1 mg) AS-21158 (5 mg)	Lot Number: See label on vial
Sequence:	H-Tyr-Pro-Tyr-Asp-Val-Pro-Asp-Tyr-Ala-OH (3-letter code) YPYDVPDYA (1-letter code)	
Molecular Weight:	1103.2	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: HA peptide is freely soluble in H₂O.

Storage: HA 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 belongs to the influenza hemagglutinin (HA) family and is responsible for attaching the virus to cell receptors and initiating infection. Ref: Schneider, C. et al. *Arch. Virol.* **125**, 103 (1992); Sleigh, MJ. et al. *J. Virol.* **37**, 845 (1981); Czech, M. et al. *J. Cell Biol.* **123**, 127 (1993).

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.*

The b_{10,11,12} HA peptide was obtained in crude form by Anaspec (San Jose, CA) and purified and characterized in the same manner. Peptides were loaded onto the purified HLA-DR1 (10–20 μM) by extended incubation (≈72 h) at 37°C using a 50- to 100-fold molar excess of peptide over HLA-DR1, in 10 mM HEPES, 50 mM NaCl, 0.02% NaN₃ (pH 7.0), in the presence of a protease inhibitor mixture consisting of 200 μM 4-(2-aminoethyl)-benzenesulfonyl fluoride, 17 μM bestatin, 2 μM E-64, 85 μM EDTA, and 20 μM pepstatin A (all from Calbiochem, San Diego, CA)- [Stratikos, E. et al. *J. Immunol.* **172**, 1109 \(2004\).](#)

Published Citations:

Stratikos, E. et al. *J. Immunol.* **172**, 1109 (2004).

Related Products:

Name	Cat#	Size
Influenza HA (307 - 319) (PKYVKQNTLKLAT)	AS-61028	1 mg
3 x Hemagglutinin (HA) Tag (MEYPYDVPDYAAEYPYDVPDYAAEYPYDVPDYAAKLE)	AS-63764	1 mg
Influenza Virus Control Peptide Pool	AS-62340	3 mg

For Research Use Only