

Product Data Sheet

Product Name: HA peptide

Catalog Number: AS-21156 (1 mg) Lot Number: See label on vial

AS-21158 (5 mg)

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 bh 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