

Product Name:	PEP1	
Catalog Number:	AS-60430-1 (1 mg)	Lot Number: See label on vial
Sequence:	Ser-Gly-Ser-Trp-Leu-Arg-Asp-Val-Trp-Asp-Trp-Ile-Cys-Thr-Val-Leu-Thr- Asp-Phe-Lys-Thr-Trp-Leu-Gln-Ser-Lys-Leu-Asp-Tyr-Lys-Asp-NH2 (3-letter code) SGSWLRDVWDWICTVLTDFKTWLQSKLDYKD-NH2 (1-letter code)	
Molecular Weight:	3806.3	

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Reconstitute by adding 100 µl 1%NH₄OH to 1 mg PEP1 peptide. Dilute this peptide solution to approximately 1 mg/ml (or more dilute) with a buffer such as PBS or another buffer; aliquot and store at -20C.

Storage: PEP1 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 synthetic peptide mimics wild-type AH (amphipathic helix) and inhibits membrane association of NS5A, hence impairing HCV replication. Ref: Glenn, JS. et al. J. Virol. 77, 6055 (2003).

Published Citations:

Chan, S. et al. Phys. Chem. Chem. Phys. 10, 3203 (2008).

Related Products:

Name

Cat # Size AS-61017-5 5 µg AS-61017-10 10 µg HCV NS3/4A protease, recombinant

For Research Use Only