



Product Data Sheet

Product Name: Hepatitis Virus C NS3 Protease Inhibitor 2

Catalog Number: AS-25346 (1 mg) **Lot Number:** See label on vial

Sequence: Ac-Asp-Glu-Dif-Glu-Cha-Cys-OH (3-letter code)
Ac-DE-Dif-E-Cha-C (1-letter code)

Molecular Weight: 914.1

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Reconstitute by adding 80-100 µl 1%NH₄OH to 1 mg Hepatitis Virus C NS3 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: Hepatitis Virus C NS3 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: A potent peptide inhibitor of Hepatitis Virus C NS3 protease. Ref: Ingallinella, P. et al. *Biochem.* **37**, 8906 (1998); Ferreon, J. et al. *J. Biol. Chem.* **280**, 20483 (2005).

Related Products:

Name	Cat #	Size
HCV (Hepatitis C Virus) NS3/4A Protease Substrate (Ac-DE-Dap(QXL520)-EE-Abu-ψ-[COO]AS-C(5-FAMsp)-NH ₂)	AS-60798	1 mg
HCV Protease FRET Substrate (RET S1) (Ac-DE-D(Edans)-EE-Abu-ψ-[COO]-AS-K(DabcyI)-NH ₂)	AS-22991 AS-22991-50 AS-22991-100	1 mg 50 mg 100 mg
HCV NS3/4A Protease Substrate Hydrolysis Product 1 (Ac-DE-Dap(QXL520)-EE-Abu)	AS-60978	1 mg

For Research Use Only