



Product Data Sheet

Product Name: HCV (Hepatitis C Virus) NS3/4A Protease Substrate

Catalog Number: AS-60798 (1 mg) **Lot Number:** See label on vial

Sequence: Ac-Asp-Glu-Dap(QXL520)-Glu-Glu-Abu-ψ-[COO]-Ala-Ser-Cys(5-FAMsp)-NH₂ (3-letter code)
Ac-DE-Dap(QXL520)-EE-Abu-ψ-[COO]AS-C(5-FAMsp)-NH₂ (1-letter code)

Molecular Weight: 1913.8

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized red powder

Peptide Reconstitution: Reconstitute by adding 100 µl 1%NH₄OH to 1 mg HCV NS3/4A Protease Substrate. Immediately dilute this peptide solution to approximately 1 mg/ml (or more dilute) with a buffer such as PBS or another buffer; aliquot into several freezer vials and store at -20C.

Storage: HCV NS3/4A 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 highly sensitive FRET substrate for assaying HCV (Hepatitis C Virus) NS3/4A protease activity. It detects < 0.1pmol of HCV NS3/4A protease. Upon cleavage of this substrate, fluorescence can be monitored at Abs/Em = 490/512 nm. Ref: Tong, XH. et al. Understanding Biology Using Peptides. Sylvie E. Blondelle(Editor): *American Peptide Society* (2005); Tang, Y. et al. *Viral Hepatitis Meeting Abstract* (2005).

Published Citations:

Seiwert, S. et al. *Antimicrob. Agents Chemother.* **52**, 4432 (2008).

For Research Use Only