

Hepatitis-C Virus

Related Products

AnaSpec, Inc. • 2149 O'Toole Ave. • San Jose, CA 95131 • 800-452-5530 • Fax 408-452-5059 • service@anaspec.com • www.anaspec.com

Product	Size	Catalog #
Peptides		
HCV Protease Substrate	250 µg	22991-025
(RET S1)Ac-DED(EDANS)EEAbu-γ-[COO]ASK(DABCYL)-NH ₂	1 mg	22991
Ex/Em=355/500 nm.	5 mg	22991-5
RET S1 Hydrolysis Product 1	1 mg	25334
Ac-DED(EDANS)-EEAbu		
RET S1 Hydrolysis Product 2	1 mg	25132-1
Lac-SK(DABCYL)-NH ₂	5 mg	25132-5
HCV (Hepatitis C Virus) NS3/4A Protease Substrate		
Ac-DE-Dap(QXL520)-EE(Abu)(COO)AS-C(5-FAMsp)-NH ₂	1 mg	60798
Ex/Em= 490/512 nm.		
HCV NS3/4A Protease Substrate Hydrolysis Product 1		
Ac-DE-Dap(QXL520)-EE(Abu)	1mg	60978
HCV NS3/4A Protease Substrate Hydrolysis Product 2		
Lac-S-C(5-FAMsp)-NH ₂	1 mg	60979
KK4A	1 mg	25335
KKGSVVIVGRIVLSGK	5 mg	24491
Pep 4A	1 mg	25336
GSVVIVGRIVLSGR-NH ₂	5 mg	24492
Pep 4AK	1 mg	25337
KKKGSVVIVGRIVLSGR-NH ₂	5 mg	24493
4A/4B Peptide (1)	1 mg	25340
DEMEECSQLPYI	5 mg	24483
4A/4B Peptide (2)	1 mg	25341
DEMEECASHLPYK(Ac)-NH ₂	5 mg	24484
4A/4B Peptide (3)	1 mg	25342
DEMEECASHLPYE(EDANS)-NH ₂	5mg	24485
4A/4B, 5A/5B Peptide	1 mg	25344
Ac-DEMEECSMSY-NH ₂	5 mg	24487
4B/5A Peptide	1 mg	25343
ECTTPCSGSWLRD	5 mg	24486
5A-pNA	1 mg	25117
EDVV-Abu-C-pNA	5 mg	25118

Product	Size	Catalog #
5A/5B Peptide (1)	1 mg	25338
EDVV-Abu-CSMSY	5 mg	24494
5A/5B Peptide (2)	1 mg	25048
Ac-EEVVAC-AMC	5 mg	25049
5A/5B Peptide (3)	1 mg	25050
Ac-EEVVAC-pNA	5 mg	25051
5A/5B Peptide (4)	1 mg	25052
K(Biotin)KEDVV-Abu-CS-Abu-SYKK-NH ₂	5 mg	25053
Ac-5A/5B Peptide	1 mg	25339
Ac-EDVVCCSMSY-NH ₂	5 mg	24482
Hepatitis Virus C NS3 Protease Inhibitor 1		
Ac-DEMEEC	1 mg	25345
	5mg	24488
Hepatitis Virus C NS3 Protease Inhibitor 2		
Ac-DEDif-E-Cha-C	1 mg	25346
	5 mg	24489
Hepatitis Virus C NS3 Protease Inhibitor 3		
Ac-DEDif-I-Cha-C	1 mg	25347
	5 mg	24490
Hepatitis Virus C NS3 Protease Inhibitor 4		
Ac-DELE-Cha-C	1 mg	25348
	5 mg	24495
HCV Substrate	1 mg	24199
EFDEMEECASHLPYI	5 mg	24411

Assay Kits

for the detection of HCV NS3/4A protease activity:

EnzoLyte™ 490 HCV Protease Assay Kit *Fluorimetric*		
(Ex/Em=355/500 nm)	1 kit	71126
EnzoLyte™ 520 HCV Protease Assay Kit *Fluorimetric*		
(Ex/Em= 490/512 nm)	1 kit	71145
EnzoLyte™ 620 HCV Protease Assay Kit *Fluorimetric*		
(Ex/Em= 591/622 nm)	1 kit	71146