

Data Sheet

HCV1b (NS3/NS4A)

Recombinant, N-terminal Flag-His tag
Catalog #: 80107
Lot #: 120817-G **Conc.:** 0.2 mg/ml

Formulated in: 50 mM Tris-HCl, pH 8.0, 500 mM NaCl, 10% glycerol, 0.2% Triton X-100, and 10 mM BME.

Stability: >6 months at -80°C

References:

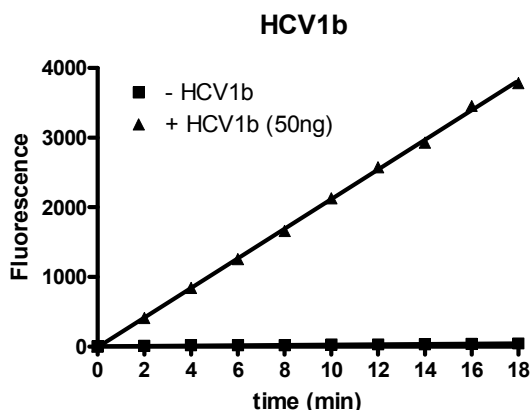
1. Lin, C., *et al.*, *Hepatitis C Viruse*. Norfolk (UK): Horizon Bioscience; 2006. Chapter 6.
2. Lang Kuhs, K.A., *et al.*, *Mol. Ther.* 2012 May; **20(3)**:669-678.
3. Massariol, M.J., *et al.*, *Biochem Biophys Res Commun.* 2010 Jan 1; **391(1)**:692-697.

Description: Fusion protein corresponding to the serine protease NS3 (a.a. 3-181) and cofactor NS4A (a.a. 21-32) from Hepatitis C virus genotype 1b (HCV1b, GenBank Accession No. AF054247), N-terminal Flag-His tag, and a 4 a.a. linker, expressed in an *E. coli* expression system. MW = 22.3 kDa.

Assay Conditions: Assay was performed using 100 µl HCV assay buffer containing 50 mM Tris, pH 7.4, 150 mM NaCl, 5 mM DTT, 10% glycerol, and 5 µM fluorescently labeled HCV substrate peptide at room temperature for 20 minutes continuously at ex350/em500.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

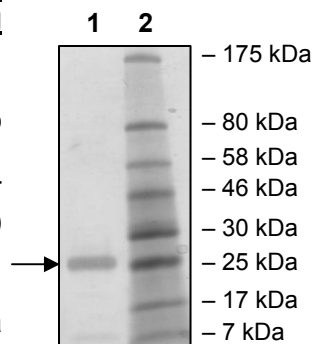
Quality Assurance



**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 2 µg HCV1b
Lane 2:
 Protein Marker
 BioLabs (#P7708L)

MW: 22.3 kDa
Purity: ≥80%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Amino Acid Sequence:

MDYKDDDDKHHHHHGSVVIVGRIILSGSGSITAYSQQTRGVLGCIITSLTGRDKNQVE
GEVQVVSTATQSFLATCINGVCWTVYHGAGSKTLAGPKGPITQMYTNVDL DLVGWQ
APPGARSMTPCSCGSSDLVTRHADVIPVRRRGDSRGSLLSPRPVSYLKGSSGGPL
LCPSGHVVGVFRAAVCTRGVAKAVDFIPVESMETTMRS

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com