

Data Sheet

HCV1a (D168V)

Recombinant, N-terminal FLAG-His tag
Catalog #: 80103
Lot #: 130808 **Conc.:** 0.084 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, and 250 mM imidazole.

Stability: >6 months at -70°C . Avoid freeze/thaw cycles.

References:

1. Lin, C., *et al.*, *Hepatitis C Virus*. Norfolk (UK): Horizon Bioscience; 2006. Chapter 6.
2. Prabhu, R., *et al.*, *Exp Mol Pathol*. 2004 Jun; **76(3)**:242-52.

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 1a (GenBank Accession No. NC_004102) with Asp-to-Val mutation on a.a. 168, and N-terminal FLAG-His tag, MW = 22.4 kDa, expressed in an *E. coli* expression system.

Application:

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

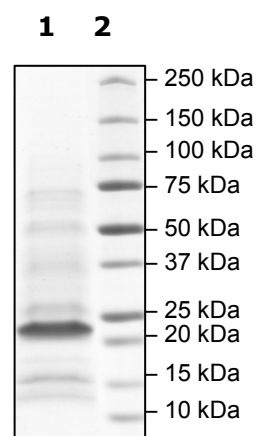
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
1.6 μg HCV1a (D168V)

Lane 2:
Protein Marker

Purity: $\geq 53\%$
MW: 22.4 kDa →



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

Sequence:

MDYKDDDDKHHHHHHGCVVIVGRIVLS**GSGS**ITAYAQQTRGLLGCIITSLTGRDKNQV
EGEVQIVSTATQTFLATCINGVCWTVYHGAGTRTIASPKGPVIQMYTNVDQDLVGWPA
PQGSRLTPCTCGSSDLVTRHADVIPVRRRGDSRGSLLSPRPISYLKGSSGGPLLC
PAGHAVGLFRAAVCTRGVAKAV**V**FIPVENLETTMRS

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