

Data Sheet

HCV1a (R155K), FLAG-His tag

Recombinant, N-terminal FLAG-His tag
Catalog #: 80101
Lot #: 130815-B **Conc.:** 0.37 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 300 mM imidazole

Stability: At least 6 months at -80°C . *Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

References:

1. Lenz, O., *et al.*, *J Hepatol.* 2013 Mar;**58(3)**:445-51.
2. Sheaffer, A.K., *et al.*, *Antivir Ther.* 2011;**16(5)**:705-18.

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 (HCV1a), GenBank Accession No. NC_004102, with Arg-to-Lys mutation on a.a. 155, 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.

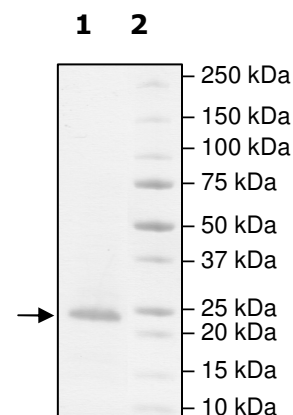
Quality Assurance

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

Lane 1:
3 μg HCV1a (R155K)

Lane 2:
Protein Marker

MW: 22.4 kDa
Purity: $\geq 90\%$



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:

MDYKDDDDKHHHHHGCVVIVGRIVLSGSGSITAYAQQTRGLLGCIITSLTGRDKNQVE
GEVQIVSTATQTFLATCINGVCWTVYHGAGTRTIASPKGPVIQMYTNVDQDLVGWPAPQ
GSRSLTPCTCGSSDLYLVTRHADVIPVRRRGDSRGSLLSPRPISYLGSSGGPLLCPAG
HAVGLFKAAVCTRGVAKAVDFIPVENLETTMRS

Red: Flag-His tag

Black without underline: NS4A cofactor (a.a. 21-32) (NP_751925)

Blue: Short linker

Black with underline: NS3 protease domain (a.a. 3-181) (NP_803144)

Yellow: R155K

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