

Data Sheet

HCV2a

Recombinant, N-terminal Flag-His tag

Catalog #: 80108

Lot #: 120927-C **Conc.:** 0.27 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 400 mM NaCl, 0.2% Triton X-100, 10 mM β -mercaptoethanol, 100 μ g/mL Flag peptide, and 30% glycerol.

Stability: >6 months at -80°C

References:

1. Jensen, T.B., *et al.*, *J. Infect. Dis.* 2008;**198**(12):1756-1765.
2. Wang, Y., *et al.*, *Cell Immunol.* 2006 May;**241**(1):38-46.

Description:

Hepatitis C virus (HCV) genotype 2a, also known as HCV2a, is a fusion protein of a.a. 21-32 of NS4A, linker (4 a.a.) and a.a. 3-181 of NS3 (GenBank Accession No. FJ393024) with N-terminal Flag-His tag, MW = 22.3 kDa, expressed in an *E. coli* expression system.

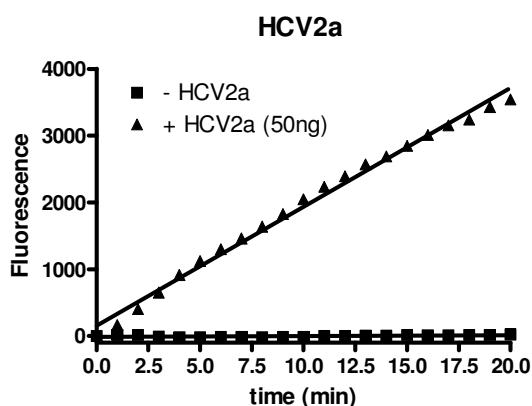
Specific Activity: 93 pmole/min/ μ g

Assay conditions: Assay was done in 100 μ l HCV assay buffer containing 50 mM Tris, pH7.4, 150 mM NaCl, 5 mM DTT, 10% glycerol, and 5 μ M HCV substrate peptide. Reaction was monitored at room temperature for 20 min continuously at ex350/em500.

Application:

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

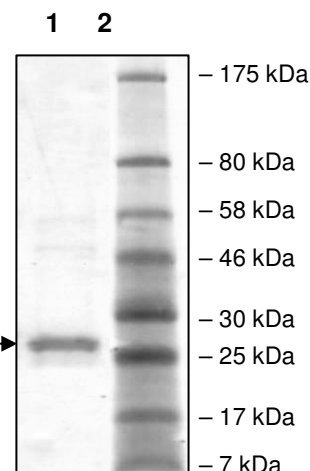
Quality Assurance



4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μ g HCV2a
Lane 2:
Protein Marker

MW: 22.3 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:

**MDYKDDDDKHHHHHHGCVSIIGRLHVNGSGSITAYAQQTRGLLGAIVVSMTGRDRTE
QAGEVQILSTVSQSFLGTTISGVLWTVYHGAGNKTLAGLRGPVTQMYSSAEGDLVGW
PSPPGTKSLEPCKCGAVDLYLVTRNADVIPARRRGDKRGALLSPRPISLTKGSSGGP
VLCPRGHVVGLFRAAVCSRGVAKSIDFIPVETLDVVTRS**

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