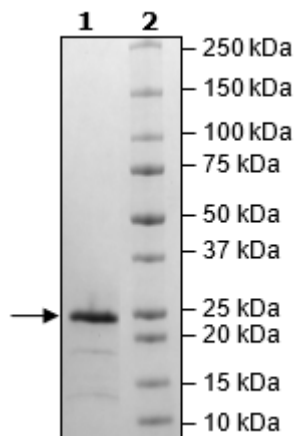


Product Information

Construct:	(FLAG-His-NS4A(21-32)-NS3(3-181)) (HCV1b)
Mutation:	NS3: R155K
Concentration:	0.28 mg/ml
Species:	Hepatitis C Virus
Formulated In:	40 mM Tris-HCl, pH 8.0, 400 mM NaCl, 0.2% Triton X-100, 10 mM beta-mercaptoethanol and 30% glycerol.
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	AF054247
MW:	22 kDa
Purity:	≥90%
Specific Activity:	≥12 pmol/min/μg
Assay Conditions:	Assay was performed in 50 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 (Anaspec #AS-22991). After 10-minute incubation at room temperature, reaction was at ex350/em500.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



HCV1b (R155K) Activity

