

**HCV NS3 genotype 1b 36 kDa**

Hepatitis C virus NS3 protein
recombinant, *E. coli*

Cat. No.	Amount
PR-1406-S	200 µg
PR-1406-L	1 mg

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Short term exposure (up to 10 days) to ambient temperature possible.

Shelf Life: 6 months

Molecular Weight: about 36 kDa

Purity: ≥ 85 %

Form: liquid (Supplied in Tris-HCl 25 mM pH 8.0, Glutathione 5 mM, Urea 1.5 M, Triton X-100 0.2 %, Glycerol 50 %)

Applications:

Suitable for use in ELISA. Each laboratory should determine an optimum working titer for use in its particular application. Other applications have not been tested but use in such assays should not necessarily be excluded.

Description:

The protein contains immunodominant region of HCV NS3, genotype 1b, amino acids 42-144. The protein is fused to a GST tag.

Specificity:

Immunoreactive with sera of HCV-infected individuals