

**HCV NS5 genotype 1b 38 kDa**

Hepatitis C virus NS5 protein  
recombinant, *E. coli*

Cat. No.	Amount
PR-1416-S	200 µg
PR-1416-L	1 mg

**For *in vitro* use only!**

**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 38 kDa

**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 NS5, genotype 1b, amino acids 61-166. The protein is fused to a GST tag.

**Specificity:**

Immunoreactive with sera of HCV-infected individuals