

**HCV NS3 genotype 2c 37 kDa**

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

Cat. No.	Amount
PR-1408-S	200 µg
PR-1408-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 37 kDa

**Purity:** ≥ 85 %

**Form:** liquid (Supplied in Sodium Carbonate-Bicarbonate 20 mM pH 10.5, Tris-HCl 50 mM, Sodium Chloride 60 mM, Glutathione 10 mM, SDS 0.1 %, 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 2c, amino acids 1361-1463. The protein is fused to a GST tag.

**Specificity:**

Immunoreactive with sera of HCV-infected individuals