

**HCV NS4 (1691-1710)(1712-1733)(1921-1940) mosaic 34 kDa**

Hepatitis C virus NS4 protein, mosaic recombinant, *E. coli*

Cat. No.	Amount
PR-1414-S	200 µg
PR-1414-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

Shelf Life: 6 months

Molecular Weight: about 34 kDa

Form: liquid (Supplied in Tris-HCl 50 mM pH 8.0, Sodium Chloride 60 mM, Glutathione 10 mM, 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 regions of HCV NS4, amino acids (1691-1710)(1921-1940)(1712-1733)(1921-1940). The protein is fused to a GST tag.

Specificity:

Immunoreactive with sera of HCV-infected individuals