

HCV-N-G5 (residues 2-119) Hepatitis C Virus Nucleocapsid Genotype 5 recombinant, *E. coli*

Cat. No.	Amount
PR-1138	100 µg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 50mM Tris-HCl pH 8.0, 60mM NaCl, 10mM glutathione, 0.25% sarkosil and 50% glycerol.

Application

Antigen in ELISA and Western blots, excellent antigen for detection of HCV with minimal specificity problems.

Specificity

Immunoreactive with sera of HCV-infected individuals.

Purity

>95% by SDS-PAGE

Description

The protein contains the HCV core nucleocapsid immunodominant regions, amino acids: 1-102.

Hepatitis C Virus core proteins are purified by proprietary chromatographic techniques.

Background

The hepatitis C virus (HCV) core protein represents the first 191 amino acids of the viral precursor polyprotein and is cotranslationally inserted into the membrane of the endoplasmic reticulum (ER)

Hepatitis C virus (HCV) core is a viral structural protein; it also participates in some cellular processes, including transcriptional regulation. However, the mechanisms of core-mediated transcriptional regulation remain poorly understood. Hepatitis C virus (HCV) core protein is thought to contribute to HCV pathogenesis through its interaction with various signal transduction pathways. In addition, HCV core antigen is a recently developed marker of hepatitis C infection. It is remarkably efficient at establishing persistent infection, suggesting that it has evolved one or more strategies aimed at evading the host immune response. T cell responses, including interferon-gamma production, are severely suppressed in chronic HCV patients. The HCV core protein has been previously shown to circulate in the bloodstream of HCV-infected patients and inhibit host immunity through an interaction with gC1qR.

Selected References:

- Kang *et al.* (2005) Proteomic profiling of cellular proteins interacting with the hepatitis C virus core protein. *Proteomics*. **5**:2227.
- Fukutomi *et al.* (2005) Hepatitis C virus core protein stimulates hepatocyte growth: Correlation with upregulation of wnt-1 expression. *Hepatology* **41**:1096.
- Gaudy *et al.* (2005) Usefulness of the hepatitis C virus core antigen assay for screening of a population undergoing routine medical checkup. *J. Clin. Microbiol.* **43**:1722.
- Lindh *et al.* (2005) Monitoring treatment response by the hepatitis C virus core antigen assay. *Eur. J. Clin. Microbiol. Infect. Dis.* **24**:230.
- Boni *et al.* (2005) Hepatitis C Virus Core Protein Acts as a trans-Modulating Factor on Internal Translation Initiation of the Viral RNA. *J. Biol. Chem.* **280**:17737.