

HAV-P3C (residues 1643-1743) Hepatitis A Virus Core Protein P3C recombinant, *E. coli*

Cat. No.	Amount
PR-1121	100 µg

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied as a 1 mg/ml solution in 10 mM CBB, pH 9.6, 0.1% SDS, and 50% glycerol.

Specificity

Immunoreactive with sera of HAV-infected individuals.

Protein synonyms/aliases

Genome polyprotein

Purity

>95% by SDS-PAGE (coomassie staining) and RP-HPLC.

Description

The protein contains the HAV probable protein P3C immunodominant regions, amino acids: 1643-1743. HAV core proteins are purified by proprietary chromatographic techniques.

Application

Recombinant HAV-P2C-P3A/2 may be used in ELISA and Western blots, excellent for detection of HAV with minimal specificity problems.

Background

Forty-two antigenic domains were identified across the hepatitis A virus (HAV) polyprotein by using a set of 237 overlapping 20-mer synthetic peptides spanning the entire HAV polyprotein and a panel of serum samples from acutely HAV-infected patients. The term "antigenic domain" is used in this study to define a protein region spanned with consecutive overlapping immunoreactive peptides.

Nineteen antigenic domains were found within the structural proteins, and 22 were found within the nonstructural proteins, with 1 domain spanning the junction of VP1 and P2A proteins. Five of these domains were considered immunodominant, as judged by both the breadth and the strength of their immunoreactivity. One domain is located within the VP2 protein at position 57-90 aa. A second domain, located at position 767-842 aa, contains the C-terminal part of the VP1 protein and the entire P2A protein. A third domain, located at position 1403-1456 aa, comprises the C-terminal part of the P2C protein and the N-terminal half of the P3A protein. The fourth domain, located at position 1500-1519 aa, includes almost the entire P3B, and the last domain, located at position 1719-1764 aa, contains the C-terminal region of the P3C protein and the N-terminal region of the P3D protein. It is interesting to note that four of the five most immunoreactive domains are derived from small HAV proteins and/or encompass protein cleavage sites separating different HAV proteins.

Selected References:

Peters *et al.* (2005) Hepatitis A virus proteinase 3C binding to viral RNA: correlation with substrate binding and enzyme dimerization. *Biochem. J.* **385**:363.

Jain R.P and Vederas J.C. (2004) Structural variations in ketoglutamines for improved inhibition against hepatitis A virus 3C proteinase. *Bioorg.*

HAV-P3C (residues 1643-1743)

Hepatitis A Virus Core Protein P3C

recombinant, *E. coli*

Med. Chem. Lett. **14**:3655.

Losick *et al.* (2003) Signals in hepatitis A virus P3 region proteins recognized by the ubiquitin-mediated proteolytic system. *Virology*. **309**:306.

Kanda *et al.* (2003) Hepatitis A virus VP3 may activate serum response element associated transcription. *Scand. J. Gastroenterol.* **38**:307.

Hu *et al.* (2002) Mutational characteristics in consecutive passage of rapidly replicating variants of hepatitis A virus strain H2 during cell culture adaptation. *World J. Gastroenterol.* **8**:872.