

Hepatitis C core protein monoclonal antibody (C7-50)

Hepatitis C virus (HCV) is the primary cause of non-A, non-B hepatitis and results in most HCV-infected people developing chronic infections, liver cirrhosis and hepatocellular carcinoma. The 21kDa core protein of HCV is well conserved among different HCV genotypes and may suppress hepatitis B virus replication in a phosphorylation dependent manner.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 10

[View Online »](#)

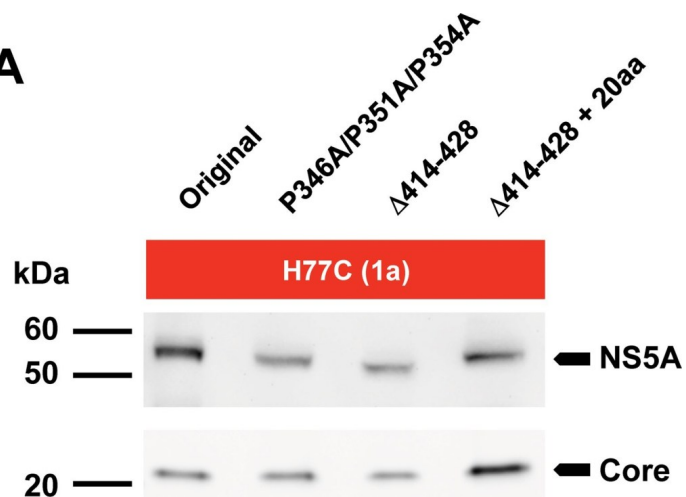
Ordering Information

[Order Online »](#)

ALX-804-277-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

A

NS5A stability of H77C(1a) mutants. Amounts of NS5A were compared to Core after spread of infection to >80% of Huh7.5 cells transfected with H77C(1a) and P346A/P351A/P354A, Δ 414-428 or " Δ 414-428+20aa" mutants. (A) Western blots from a representative experiment of NS5A (56/58 kDa for original), and Core (21 kDa) detected with antibodies H26 (Abcam) and C7-50 (Enzo Life Science), respectively. (B) Ratios between NS5A/Core were calculated after subtraction of background using ImageJ and normalized to the original H77C(1a). Values from two independent transfection experiments are shown.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Analysis of functional differences between hepatitis C virus NS5A of genotypes 1-7 in infectious cell culture systems. *PLoS Pathog* (2012)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA, ICC, IP, WB

Application Notes Immunocytochemical staining of HCV core protein in transfected chimp hepatocytes yields a staining pattern consistent with cytoplasmic and vesicular staining. Detects a band of ~21kDa by Western blot.

Clone C7-50

Formulation Liquid. In PBS containing 0.05% sodium azide.

Host Mouse

Immunogen Purified HCV (hepatitis C virus) core fusion protein (genotype 1b).

Isotype IgG1

Recommendation Dilutions/Conditions Western Blot (1µg/ml) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Hepatitis C virus

Specificity Recognizes an epitope between aa 21-40 of HCV core protein. This sequence is conserved among different HCV strains.

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)