

HAV-VP1-P2A

(residues 669-782)

Ordering Information

[Order Online »](#)

JBS-PR-1114	100µg
-------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Hepatitis A Virus Coat Protein VP1-Core Protein P2A

Application Notes Recombinant rHAV-VP1-P2A/Mosaic Antigen may be used in ELISA and Western blots, excellent for detection of HAV with minimal specificity problems.

Formulation liquid (Supplied in 10 mM Tris-HCl pH 9.6, 1.5 M urea and 50% glycerol)

MW 23.5 kDa

Purity > 90 % (SDS-PAGE)

Source recombinant, E. coli

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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