

Hepatitis B Virus

Surface Antigen – adr subtype

Ordering Information

[Order Online »](#)

JBS-PR-1197	100µg
-------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling do not freeze

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes Antigen in ELISA, Immunochromatography and monoclonal antibody production.

Formulation liquid (Supplied in 20 mM sodium phosphate buffer (PB), 154 mM NaCl, pH 7.1)

MW 23 – 27 kDa

Purity > 95 % (RP-HPLC)

Source recombinant, Chinese hamster ovary (CHO) cells

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-be@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