

West Nile Virus Pre-M His

Ordering Information

[Order Online »](#)

JBS-PR-1274	100µg
-------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	WNV Pre-M
Application Notes	Antigen in ELISA and Western Blots, excellent antigen for detection of West-Nile virus with minimal specificity problems.
Formulation	liquid (Supplied in 20 mM Phosphate buffer pH 7.5)
Purity	> 95 % (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-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