

TmpA partial (23-41aa and 288-325aa)

Ordering Information

[Order Online »](#)

JBS-PR-1269

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 Treponema Pallidum Membrane Protein A

Application Notes Treponema Pallidum protein binds to murine anti-TmpA protein monoclonal antibodies and Treponema pallidum converted human serum polyclonal antibodies in ELISA and Western Elisa. For Dot Blots and Lateral Flow immunochromatographic diagnostic tests.

Formulation liquid (Supplied in 10 mM Tris-HCl pH 8.0, 20 mM DDT, 1 mM EDTA and 8 M urea)

MW 68 kDa

Purity > 90 % (SDS-PAGE)

Source Treponema pallidum, recombinant, E. coli

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

image not found or type not known
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