

CMV pp65 (residues 297-510)

Ordering Information

[Order Online »](#)

JBS-PR-1251	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 Cytomegalo Virus Phosphoprotein 65

Application Notes Antigen in ELISA and Western blots, excellent antigen for detection of CMV with minimal specificity problems.

Formulation liquid (Supplied in 50 mM Tris-HCl, 10 mM Glutathione, 60 mM NaCl, 0.125 % Sarcosil and 50% glycerol)

MW 52.2 kDa +26 kDa GST

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