

Lom-TK 2 polyclonal antibody

Ordering Information

[Order Online »](#)

JBS-ABD-045	100µl
-------------	-------

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 Locusta-Tachykinin 2

Application ELISA, ICC

Formulation liquid (Supplied as serum, preserved in glycerol) This antiserum is obtained from rabbits after immunization with the conjugates: Loc-Tachykinin 2-glutardialdehyde-thyroglobulin.

Host Rabbit

Immunogen Synthetic peptide, APLSGFYGVR-NH₂, corresponding to the whole insect neuropeptide Locusta-Tachykinin 2 (Lom-TK 2).

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