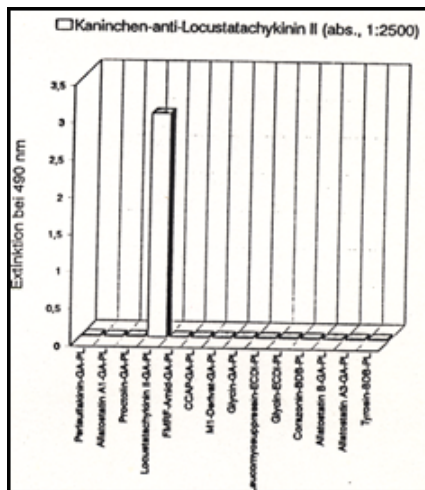




anti-Lom-TK 1

anti-Locusta-Tachykinin 1
rabbit, polyclonal

Cat. No.	Amount
ABD-044	100 µl



The two isoforms LomTK 1 and 2 have identical acarboxy-terminus heptapeptides. The anti-LomTK 2 finds out the LomTK 1 antigen.

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Form: liquid (Supplied as serum, preserved in glycerol (50 % saturation))

This antiserum is obtained from rabbits after immunization with the conjugates: Loc-Tachykinin 2-glutaraldehyde-BSA.

Applications:

Anti-Lom-TK 1 can be used for ELISA and Immunocytochemistry.

Description:

Polyclonal antibody to Locusta-Tachykinin 1.

Immunogen:

Synthetic peptide, APLSGFYGVR-NH₂, corresponding to the whole insect neuropeptide Locusta-Tachykinin 1 (LomTK 1).

Specificity:

Rabbit anti-LomTK 1 specificity is performed by an ELISA using different insect neuropeptide conjugates: Perisulfakinin-G-PL, Allatostatin A1-G-PL, Allatostatin A3-G-PL, Proctolin-G-PL, Locusta-tachykinin 2-G-PL

FMRF-amid-G-PL, CCAP-G-PL, M1-G-PL, Leucomyosuppressin-G-PL, Corazonin-G-PL

(G= glutaraldehyde, PL = polylysine).

Selected References:

Clynen *et al.* (2009) Peptidomic survey of the locust neuroendocrine system. *Insect Biochem Mol Biol.* 39:491.

Neupert *et al.* (2009) Neuropeptides in Heteroptera: identification of allatotropin-related peptide and tachykinin-related peptides using MALDI-TOF mass spectrometry. *Peptides* 30:483.

Veenstra *et al.* (1995) Immunohistological localization of regulatory peptides in the midgut of the female mosquito *Aedes aegypti*. *Histochem Cell Biol.* 104:337.

Agricola *et al.* (1995) Comparative aspects of petidergic signaling pathways in the nervous systems of arthropods. In the *Nervous system of Invertebrates*. Ed. Breidbach & Kutsch, BirkhäuserVerlag. Chung *et al.* (2010) The *Drosophila* homolog of methionine sulfoxide reductase A extends lifespan and increases nuclear localization of FOXO. *FEBS Lett.* 584: 3609.