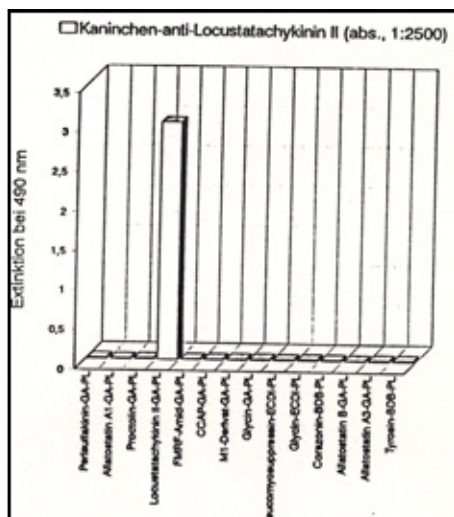


anti-Lom-TK 2 anti- Locusta-Tachykinin 2 rabbit, polyclonal antibody

Cat. No.	Amount
ABD-045	100 µl

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles



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

Form

Liquid. This antiserum is obtained from rabbits after immunization with the conjugates: Loc-Tachykinin 2-glutarialdehyde-thyroglobulin. Supplied as serum, preserved in glycerol (50% saturation).

Application

Anti- Lom-TK 2 can be used for ELISA and Immunocytochemistry.

Description

Polyclonal antibody to Locusta-Tachykinin 2

Immunogen

Synthetic peptide, APLSGFYGVN-NH₂, corresponding to the whole insect neuropeptide Locusta-Tachykinin 2 (LomTK 2)

Specificity

Rabbit anti-LomTK 2 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= glutardialdehyde, 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 peptidergic 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.