



anti-LMS

anti-Leucomyosuppressin
rabbit, polyclonal

Cat. No.	Amount
ABD-059	100 µl

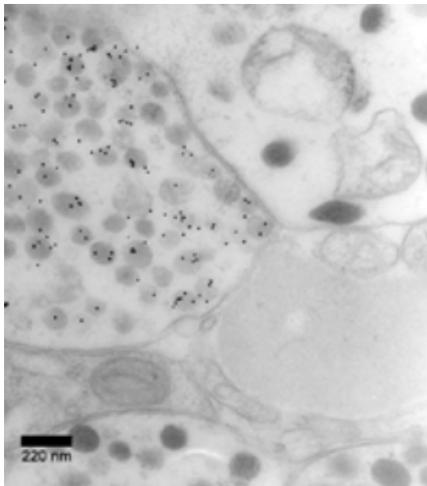


Fig. 1: The distribution of LMS-immunoreactivity in the Corpus cardiacum of male adult *Periplaneta americana*. Postembedding immunogold labelling for LMS revealed that the majority of the gold particle was bound to one type of neurosecretory granules.

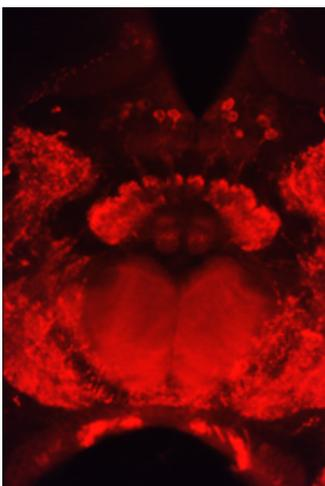


Fig. 2

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)

Applications:

Anti Leucomyosuppressin can be used for ELISA and Immunocytochemistry.

Description:

Leucomyosuppressin, a member of the FMRF amide family first characterized from the cockroach *L. maderae*. This peptide has subsequently been identified in other insects, suggesting that it is highly conserved. Leucomyosuppressin (LMS) is an inhibitory modulator of heartbeat frequency in arthropods. The anti Leucomyosuppressin serum was raised in rabbits against the conjugate from synthetic pEDVDHVFLRFamide-glutaraldehyde-thyroglobuline.

Specificity:

The anti Leucomyosuppressin serum displays no cross-reactivity with the insect neuropeptides Allatostatin, Corazonin, CCAP, Locustatachykinin II and Proctolin as tested by non-competitive ELISA. A crossreactivity to other myosuppressins (X1DVX2HX3FLRFamide) is untried. Postembedding immunogold labelling for LMS and Perisulfakinin (other RFa like peptide) revealed the gold particle were bound to different types of neurosecretory granules (Fig. 1). Immunocytochemical methods of brain (Fig. 2) and ventral nerve cords of *Periplaneta americana* indicate a specific distribution pattern.

Selected References:

Maestro *et al.* (2011) Leucomyosuppressin modulates cardiac rhythm in the cockroach *Blattella germanica*. *J Insect Physiol.* 57:1677.

Holman *et al.* (1986) Isolation, primary structure and synthesis of leucomyosuppressin, an insect neuropeptide that inhibits spontaneous contractions of the cockroach hindgut. *Comparative Biochemistry and Physiology - C Pharmacology Toxicology and Endocrinology* 85:329.