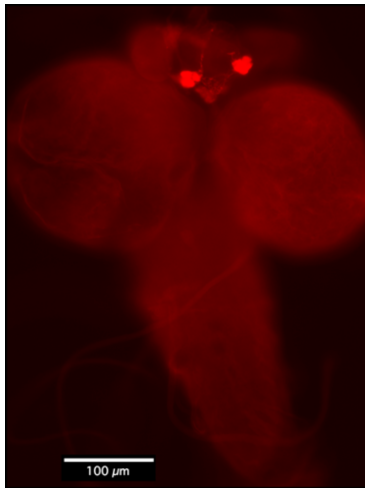


**anti-AKH I**

anti-Adipokinetic Hormone I
rabbit, polyclonal

Cat. No.	Amount
ABD-063	100 µl



Ventral nerve cord of the third larval stage from *Drosophila melanogaster*. Note the immunoreactive cells (red spots) exclusively in endocrine cells of the corpora cardiaca (ring gland).

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 AKH I can be used for ELISA and Immunocytochemistry.

Description:

Peptides belonging to the AKH/ RPCH family are produced in cells of either the corpora cardiaca (CC) of insects or of the X-organ cells in the eyestalks of decapod crustaceans. The adipokinetic hormone I (AKH I) (also named neuro-hormone D, Baumann and Penzlin, 1984) regulate energy homeostasis in insects. The hormone mobilizes energy reserves from storage, particular from fat body. There is enough evidence to show that AKH regulates the balance of carbohydrate and lipid oxidation in flight muscles.

The anti-adipokinetic hormone I serum was raised in rabbits against the conjugate from synthetic PVNFSPNWamide glutaraldehyde-thyroglobuline.

Specificity:

The anti-AKH I serum displays no cross-reactivity with insect neuropeptides Allatostatin, Corazonin, CCAP, FMR Famide, Leu-comyosuppressin, Locusta-Tachykinin 2, Perisulfakinin, and Proctolin as tested by non-competitive ELISA. Immunocytochemical methods of the ventral nerve cords of *Periplaneta americana* revealed a specific distribution of AKH-immuno-reactivity exclusively in endocrine cells of the corpus cardiacum but not seen in brain or other ganglia. In conclusion, AKHs are not synthesized in ventral nerve cords of insects.

Selected References:

Wicher *et al.* (2006) Differential receptor activation by cockroach adipokinetic hormones produces differential effects on ion currents, neuronal activity, and locomotion. *J Neurophysiol.* 96:965.

Van der Horst *et al.* (2001). Adipokinetic hormones of insect release, signal transduction, and responses. *Int Rev Cytol.* 211: 179.

Baumann *et al.* (1984) Sequence analysis of neurohormone D, a neuropeptide of an insect, *Periplaneta americana*. *Biomed Biochim Acta.* 43:K13.