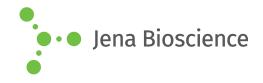
DATA SHEET





anti-PVK anti-Periviscerokinin rabbit, polyclonal

Cat. No.	Amount
ABD-065	100 μl

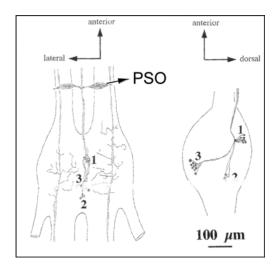


Fig. 1



Fig. 2: Example for the specificity of the anti-periviscerokinin-serum. Note the immunoreactive neurons in the ventral nerve cord of larval *Drosophila melanogaster*.

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-Periviscerokinin can be used for ELISA and Immunocytochemistry.

Description:

The anti-periviscerokinin serum was raised against the Periplaneta americana periviscerokinin-1 (Predel et.al. 1995), GASGLIPVMRNamide (PVK-1), coupled to bovine thyroglobulin using glutaraldehyde. PVK-1 is one of 24 different neuropeptides of the FMR Famide gene (Predel et.al. 2004).

Specificity:

In agreement with the mass spectrometric methods (Predel et al.2004) revealed immunocytochemical methods a tagma specific distribution of periviscerokinin-1-immunoreactivity exclusively in the abdominal perisympathetic organs but never in thoracic PSO or the retrocerebral complex of Periplaneta americana (Fig 1: PSO of third abdominal ganglia). In the contrary a anti-PVK-1-immunoreactivity is displayed in the thoracic PSOs of the ventral nerve cord from larval Drosophila melanogaster (Fig 2).

Selected References:

Predel *et al.* (2004) Unique accumulation of neuropeptides in an insect: FMRFamide-related peptides in the cockroach, Periplaneta americana. Eur J Neurosci. 6:1499.

Predel *et al.* (1995) Periviscerokinin (Pea-PVK): A novel myotropic neuropeptide from the perisympathetic organs of the American cockroach. Peptides 16: 61.