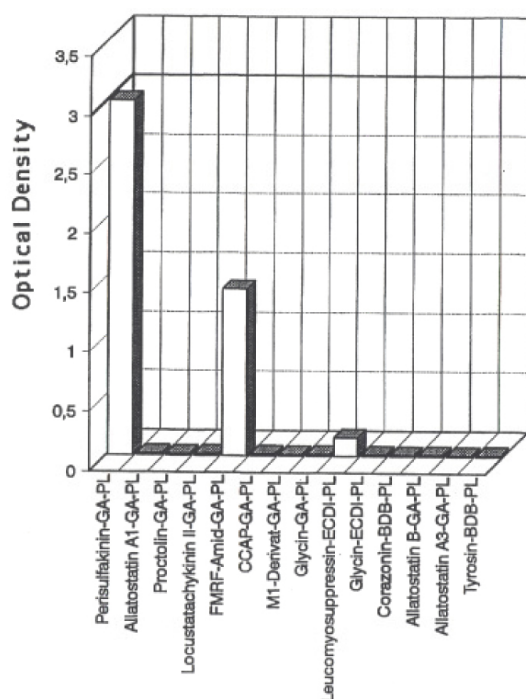


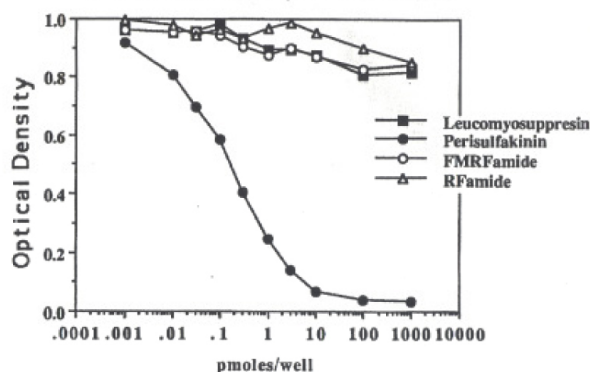
anti-PSK
anti-Perisulfakinin
rabbit, polyclonal

Cat. No.	Amount
ABD-037	100 μ l

□ rabbit-anti-perisulfakinin (1 : 20.000)



□ rabbit-anti-perisulfakinin



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

Avoid freeze / thaw cycles

Form

Liquid. Supplied as serum and 50% glycerol. These antisera are obtained from rabbits after immunization with Perisulfakinin-Glutaraldehyd-Polylysine conjugates.

Description

Polyclonal antibody to insect Perisulfakinin.

Immunogen

Synthetic peptide: EQFDDY(SO₃)GHMRF-Amid (Bachem AG/Switzerland) corresponding to the whole insect neuropeptide Perisulfakinin.

Specificity

Rabbit anti-Perisulfakinin specificity was performed by an ELISA using different insect neuropeptide conjugates: Perisulfakinin-G-PL, Allatostatin A1-G-PL, Allatostatin A3-G-PL, Proctolin-G-PL, LocustatachykininII-G-PL, FMRF-amid-GA-PL, CCAP-G-PL, M1-G-PL, Leucomyosuppressin-G-PL, Corazonin-G-PL. A weekly cross-reactivity to other RF-neuropeptides are visible (see ELISA test)

Application

Immunocytochemistry and ELISA

Selected References:

Wicher *et al.* (2007) The satiety signaling neuropeptide perisulfakinin inhibits the activity of central neurons promoting general activity. *Front Cell Neurosci.* **1**:3 Epub.
 Agricola *et al.* (1995) Comparative aspects of peptidergic signaling pathways in the nervous systems of arthropods. *EXS* **72**:303.