





Cat No. Amount

# anti-PSKR

anti-Perisulfakinin Receptor rabbit, polyclonal

Cat. No.	Amount
ABD-038	100 µl
	A thoracic M ganglia abdominal ganglia M pSKR - 434bp - 267bp
	B M DUM neurons M - 434bp - 267bp
	- 434bp - 267bp
	→ 67.8 kDa 50 kDa NS G
	-

# 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 and glycerol) These antisera are obtained from rabbits after immunization with AIKETERHKCR peptide motif (Biotrend, Colgne, Germany).

## **Applications:**

Immunocytochemistry, ELISA and Western Blots.

# Description:

Polyclonal antibody to insect Perisulfakinin Receptor.

### Specificity:

RT-PCR and Western blot analysis of PSKR expression. (A) Nested RT-PCR analysis of the PSKR expression in brain and ganglia. A 362 bp PCR product was amplified in all neuronal samples tested. (B) Single-cell RT-PCR analysis of isolated DUM neurons. In 7 out of 10 DUM neurons a PSKR-specific band was found (upper panel). As an endogenous expression control an actin-specific RT-PCR was used (lower panel). M, marker pBR322 DNA BsuRI digested (two marker bands are indicated). (C) Western blot analysis of PSKR expression in nervous system (NS) and gut (G). A 68 kDa band was found in gut only.

#### Immunogen:

Synthetic peptide motif AIKETERHKCR (Biotrend, Colgne, Germany)located in the loop between TMS4 and TMS5.

#### 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.

