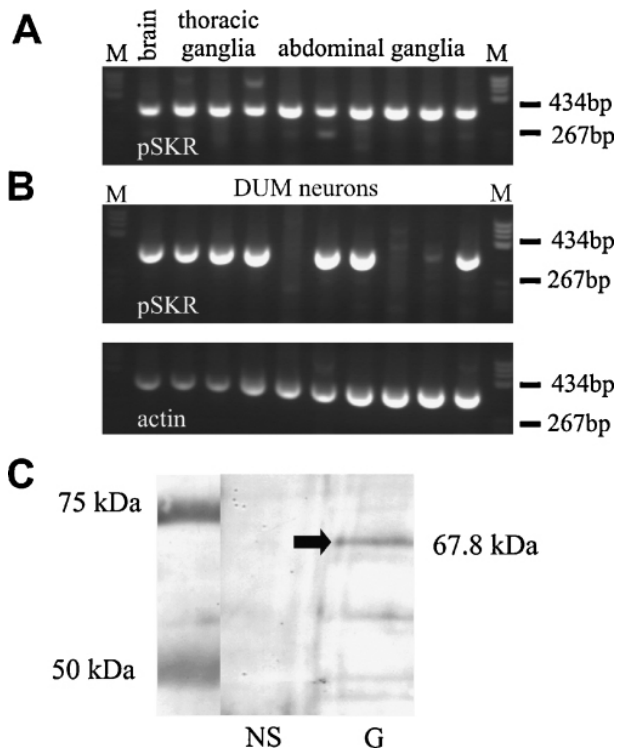


## anti-PSKR anti-Perisulfakinin Receptor rabbit, polyclonal

Cat. No.	Amount
ABD-038	100 $\mu$ l



### Description

Polyclonal antibody to insect Perisulfakinin Receptor.

### Immunogen

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

### 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.

### Application

Immunocytochemistry, ELISA and Western Blots.

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 AIKETERHKCR peptide motif (Biotrend, Colgne, Germany).

### 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.