# **SNAP 25 (CT)**polyclonal antibody

SNAP-25 is a SNAP receptor (SNARE) protein which complexes with syntaxin and synaptobrevin-2 (VAMP) to mediate membrane docking and fusion.

This antibody is covered by our Worry-Free Guarantee.

## **Ordering Information**

**Order Online** »

BML-SL3730-0025

25µl

Manuals, SDS & CofA

View Online »

### **Handling & Storage**

Long Term Storage -20°C

Shipping Blue Ice

# Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name Synaptosome-associated protein of 25kDa

**Application** ELISA, IHC, WB

**Formulation** Liquid. Neat serum containing 0.1% sodium azide.

GenBank ID M22012

**Host** Rabbit

**Immunogen** Synthetic heptadecapeptide corresponding to aa

C190RIDEANQRATKMLGSG206 of the C-terminal of

mouse SNAP-25.

Species Reactivity Bovine, Rat

UniProt ID P60879

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024