

# 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](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)