Somatostatin polyclonal antibody

Since the discovery of somatostatin-14 (SOM-14) in ovine hypothalamus, several other forms of somatostatin-like immunoreactants of different molecular weight have been reported, including somatostatin-28 which expresses the full SOM-14 sequence at its C-terminus (SOM-28[15-28]). In mammals both molecular forms are encoded at the C-terminus of the somatostatin precursor molecule. Hundreds of immunolocalisation studies have been published demonstrating the existence of somatostatin-like molecules in neurons, in pancreatic islet D cells and in endocrine cells distributed throughout the mammalian gut, as well as in rare endocrine tumours (so-called somatostatinomas). The molecular forms appear to be co-packaged in dense-cored secretory granules, frequently with other transmitter substances. A recent paper has shown that the majority of somatostatin-containing axons in the rat spinal cord express vesicular gluatamate transporter protein-2 (VGLUT2).

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

Ordering Information

Order Online »

BML-SZ1116-0025

25µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Dilute to working strength with 50mM PBS (pH 7.4) containing 1.5% sodium chloride

and 1% normal goat serum (if a goat anti-rabbit IgG linker antibody is to be used). Store

diluted antibody at 2-4°C (do not freeze) and use within 1 month.

Handling Avoid freeze/thaw cycles. Aliquot undiluted antibody into smaller volumes (not less than

10µI) prior to freezing if appropriate. The use of high quality 'antiserum-grade' plastic or

glass vials is recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC (FS), IHC (PS)

Formulation Liquid serum which has been clarified, but not purified, and

contains 0.01M sodium azide.

Host Rat

Immunogen Somatostatin-28.

Species Reactivity Species independent

Specificity Recognizes somatostatin in a wide range of species.

UniProt ID P61278

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

.

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com