

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 glutamate 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 10µl) 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 .



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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