

Somatostatin 14 polyclonal antibody

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

BML-SA1267-0025	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 SOM-14

Application IHC, Immunoelectron microscopy

Crossreactivity Does not cross-react with the N-terminus of somatostatin-28.

Formulation Liquid. Neat serum containing 0.01M sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 15-28 of human somatostatin-14.

Species Reactivity Species independent

Specificity Recognizes the C-terminus of somatostatin-14 in a wide range of mammalian and non-mammalian species, including mouse and human.

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