

Serotonin polyclonal antibody

Serotonergic cell bodies have been identified in the CNS of many vertebrate species, their processes being found widely spread throughout the brain and the spinal cord. Serotonin (5-HT)-containing fibres are found innervating the gut in many vertebrates and 5-HT is also co-stored with biologically active peptides in elements of the neuroendocrine system. Enterochromaffin (EC) cells contain high levels of serotonin and this cell type is found throughout the vertebrate gut mucosa.

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

Ordering Information

[Order Online »](#)

BML-SA1264-0025	25µl
BML-SA1264-0100	100µl

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 5-Hydroxytryptamine, 5-HT

Application IHC

Formulation Liquid. Neat serum containing 0.1% sodium azide.

Host Rabbit

Immunogen Serotonin-glutaraldehyde-BSA conjugate.

Species Reactivity Species independent

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)