## Serotonin polyclonal antibody

Serotoninergic 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	25μΙ
BML-SA1264-0100	100μΙ

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

**IHC Application** 

**Formulation** Liquid. Neat serum containing 0.1% sodium azide.

Rabbit Host

**Immunogen** Serotonin-glutaraldehyde-BSA conjugate.

**Species Reactivity** Species independent

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

Last modified: May 29, 2024

