

Serotonin Goat polyclonal antibody

Ordering Information

[Order Online »](#)

IMS-IS1035	50µl
------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Store at +4°C for short term (1-2 months). Aliquot and store at -20°C for long term. Avoid repeated freeze / thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	5-HT, 5-hydroxytryptamine
Application	IF, IHC
Application Notes	Optimized protocol with STAINperfect immunostaining kit A
Formulation	Liquid. In PBS containing 0.09% sodium azide.
Host	Goat
Immunogen	Conjugated 5-HT
Recommendation Dilutions/Conditions	1/250 -1/1000 dilution – STAINperfect
Species Reactivity	Species independent
Technical Info / Product Notes	For the Original Manufacturer's data sheet please click here .



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