

Anti-conjugated Serotonin antibody

Ref: IS1007

Clonality	Polyclonal antibody
Host	Rabbit
Applications	<u>IHC</u> / IF
Species reactivity	Reacts with all species
References	Not yet cited to our knowledge.
Format	50µl

Product information

Product overview

Product name	Anti-conjugated Serotonin rabbit polyclonal antibody
Synonyms	Anti-5-HT rabbit polyclonal antibody Anti-5-hydroxytryptamine rabbit polyclonal antibody
Immunogen	Conjugated 5-HT
Specificity	When tested in competitive ELISA, the anti-conjugated Serotonin antibody did not show any significant cross reactivity with 5-Hydroxytryptamine analogs, including Tryptamine and 5-Methoxytryptamine.
Lot number	140201

Reconstitution & storage

Form	Lyophilized powder
Purity	Purified anti-serum
Storage	Store at 4 °C
Storage buffer	Before use, vial should be resuspended in 50 µL of ultrapure water. Store at +4 °C for short term (1-2 weeks). Aliquot and store at -20 °C for long term. Avoid repeated freeze / thaw cycles.

Applications

Immunohistochemistry (IHC)	Optimal working conditions must be determined by the end-user.
-----------------------------------	--

Immunofluorescence (IF)	Optimal working conditions must be determined by the end-user.
--------------------------------	--

Restrictions	For research use only
---------------------	-----------------------

References

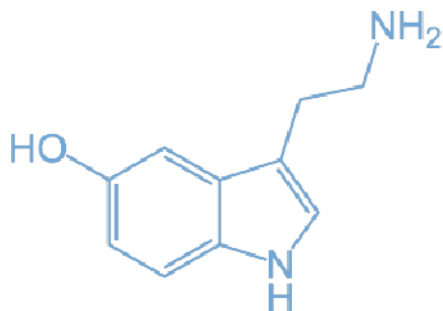
Antibody not yet cited.

Selected papers about serotonin:

- [Gershon MD. 5-Hydroxytryptamine \(serotonin\) in the gastrointestinal tract. Curr Opin Endocrinol Diabetes Obes. 2013 Feb;20\(1\):14-21. doi: 10.1097/MED.0b013e32835bc703.](#)
- [Baganz NL, Blakely RD. A dialogue between the immune system and brain, spoken in the language of serotonin. ACS Chem Neurosci. 2013 Jan 16;4\(1\):48-63. doi: 10.1021/cn300186b. Epub 2012 Dec 7.](#)
- [Bortolato M et al. The role of the serotonergic system at the interface of aggression and suicide. Neuroscience. 2013 Apr 16;236:160-85.](#)
- [Kiyasova V et al. A genetically defined morphologically and functionally unique subset of 5-HT neurons in the mouse raphe nuclei. J Neurosci. 2011 Feb 23;31\(8\):2756-68. doi: 10.1523/JNEUROSCI.4080-10.2011.](#)

Pictures

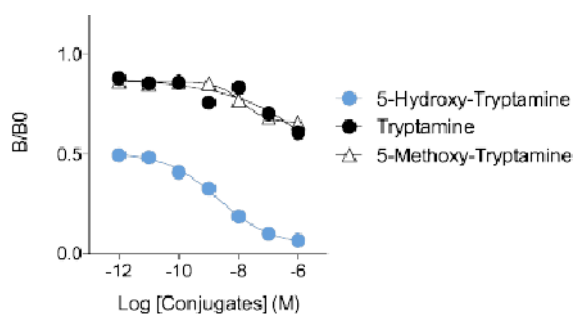
Serotonin - 5-hydroxytryptamine (5-HT)



Serotonin

Affinity & specificity of anti-5-HT antibody

Competitive ELISA demonstrates that low amounts of Serotonin conjugate are required to abolish antigen-antibody reaction (high affinity), while rising concentrations of Tryptamine and 5-Methoxy-Tryptamine conjugates do not affect reaction (high specificity).



distributed in the US/Canada by:

EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

www.EagleBio.com • info@eaglebio.com



For further information about this kit, its application or the procedures in this kit insert, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.