

Anti-conjugated Dopamine antibody

Ref: IS1005

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

Product information

Product overview

Product name	Anti-conjugated Dopamine antibody - Rabbit pAb
Synonyms	Anti-Dopamine polyclonal antibody Anti-3,4-dihydroxyphenethylamine polyclonal antibody
Immunogen	Conjugated Dopamine
Specificity	When tested in competitive ELISA, the anti-conjugated Dopamine antibody did not show any significant cross reactivity with Tyramine and L-Dopa conjugates.
Lot number	140301

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

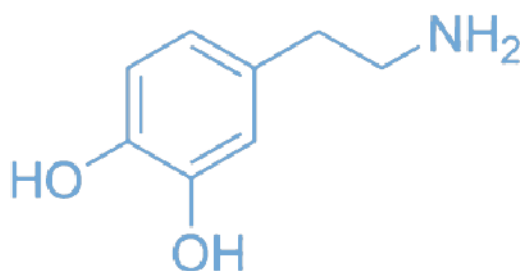
References

Antibody not yet cited.

- [Espinosa-Oliva AM et al. Role of dopamine in the recruitment of immune cells to the nigro-striatal dopaminergic structures. Neurotoxicology. 2014 Mar;41:89-101. doi: 10.1016/j.neuro.2014.01.006. Epub 2014 Jan 30.](#)
- [Watabe-Uchida M, Zhu L, Ogawa SK, Vamanrao A, Uchida N. Whole-brain mapping of direct inputs to midbrain dopamine neurons. Neuron. 2012 Jun 7;74\(5\):858-73. doi: 10.1016/j.neuron.2012.03.017.](#)
- [Chaudhury D et al. Rapid regulation of depression-related behaviours by control of midbrain dopamine neurons. Nature. 2013 Jan 24;493\(7433\):532-6. doi: 10.1038/nature11713. Epub 2012 Dec 12.](#)
- [Salamone JD, Correa M. The mysterious motivational functions of mesolimbic dopamine. Neuron. 2012 Nov 8;76\(3\):470-85. doi: 10.1016/j.neuron.2012.10.021.](#)
- [Threlfell S, Lalic T, Platt NJ, Jennings KA, Deisseroth K, Cragg SJ. Striatal dopamine release is triggered by synchronized activity in cholinergic interneurons. Neuron. 2012 Jul 12;75\(1\):58-64. doi: 10.1016/j.neuron.2012.04.038.](#)
- [Tritsch NX, Ding JB, Sabatini BL. Dopaminergic neurons inhibit striatal output through non-canonical release of GABA. Nature. 2012 Oct 11;490\(7419\):262-6. doi: 10.1038/nature11466. Epub 2012 Oct 3.](#)

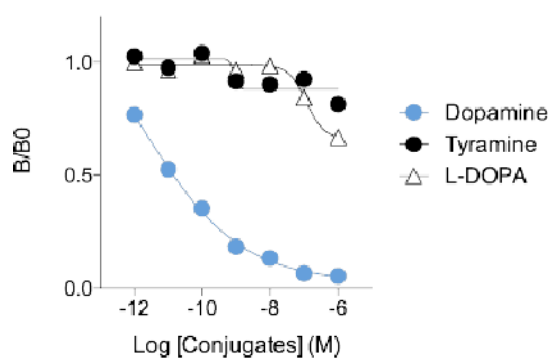
Pictures

Dopamine (3,4-dihydroxyphenethylamine)



Dopamine

Anti-Dopamine pAb (Rabbit)



Affinity & specificity of anti-Dopamine antibody

Competitive ELISA demonstrates that low amounts of Dopamine conjugate are required to abolish antigen-antibody reaction (high affinity), while rising concentrations of Tyramine and L-Dopa 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



EAGLE
BIOSCIENCES

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.