



Anti-conjugated Xanthurenic acid pAb

Ref: IS1014

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

Product information

Product overview

Product name	Anri-conjugated Xanthurenic acid rabbit polyclonal antibody
Synonyms	Anti-Xanthurenic acid rabbit polyclonal antibody 8-Hydroxy-4-oxo-1H-quinoline-2-carboxylic acid rabbit polyclonal antibody Xanthuric acid rabbit polyclonal antibody Xanthurenate rabbit polyclonal antibody 8-Hydroxykynurenic acid rabbit polyclonal antibody
Immunogen	Conjugated Xanthurenic acid
Specificity	When tested in competitive ELISA, the anti-conjugated Xanthurenic acid antibody did not display any significant cross-reactivity with analog Kynurenic acid conjugate.
Lot number	140801

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

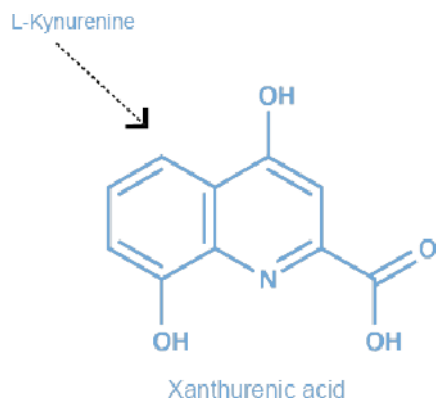
Antibody not yet cited.

Selected publications on Xanthurenic acid:

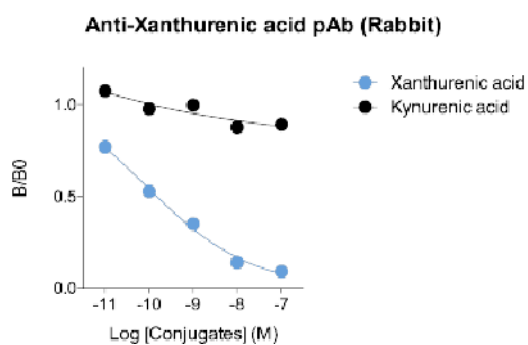
- [Neale SA, Copeland CS, Salt TE. Effect of VGLUT inhibitors on glutamatergic synaptic transmission in the rodent hippocampus and prefrontal cortex. *Neurochem Int.* 2013 Oct 10. pii: S0197-0186\(13\)00250-7. doi: 10.1016/j.neuint.2013.10.001. \[Epub ahead of print\]](#)
- [Neale SA et al. Modulation of hippocampal synaptic transmission by the kynurenine pathway member xanthurenic acid and other VGLUT inhibitors. *Neuropsychopharmacology.* 2013 May;38\(6\):1060-7. doi: 10.1038/npp.2013.4. Epub 2013 Jan 7.](#)
- [Copeland CS1, Neale SA, Salt TE. Actions of Xanthurenic acid, a putative endogenous Group II metabotropic glutamate receptor agonist, on sensory transmission in the thalamus. *Neuropharmacology.* 2013 Mar;66:133-42. doi: 10.1016/j.neuropharm.2012.03.009. Epub 2012 Apr 2.](#)
- [Taleb O et al. Xanthurenic acid binds to neuronal G-protein-coupled receptors that secondarily activate cationic channels in the cell line NCB-20. *PLoS One.* 2012;7\(11\):e48553. doi: 10.1371/journal.pone.0048553. Epub 2012 Nov 6.](#)

Pictures

Xanthurenic acid (XA)



Affinity & specificity of Xanthurenic acid polyclonal antibody



Competitive ELISA demonstrates that low amounts of Xanthurenic acid conjugate are required to abolish antigen-antibody reaction (high affinity), while rising concentrations of Kynurenic acid conjugate 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.