



Anti-conjugated picolinic acid antibody

Ref: IS1015

Clonality	Polyclonal
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 Picolinic acid rabbit polyclonal antibody
Synonyms	pyridine-2-carboxylic acid polyclonal antibody
Immunogen	Conjugated picolinic acid
Specificity	When tested in competitive ELISA, the anti-Picolinic acid antibody did not show any significant cross reactivity with analog Quinolinic acid.
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

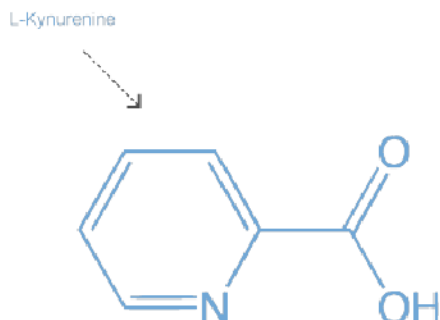
Antibody not yet cited.

Selected publications about Picolinic acid:

- [Zavitsanou K et al. Effect of maternal immune activation on the kynurenine pathway in preadolescent rat offspring and on MK801-induced hyperlocomotion in adulthood: Amelioration by COX-2 inhibition. Brain Behav Immun. 2014 May 27. pii: S0889-1591\(14\)00133-0. doi: 10.1016/j.bbi.2014.05.011.](#)
- [Reyes Ocampo J et al. Kynurenines with neuroactive and redox properties: relevance to aging and brain diseases. Oxid Med Cell Longev. 2014;2014:646909. doi: 10.1155/2014/646909. Epub 2014 Feb 17.](#)
- [Zuwala-Jagiello, J et al. Picolinic Acid in Patients with Chronic Hepatitis C Infection: A Preliminary Report. Mediators Inflamm. 2012;2012:762863. doi: 10.1155/2012/762863. Epub 2012 Apr 10.](#)

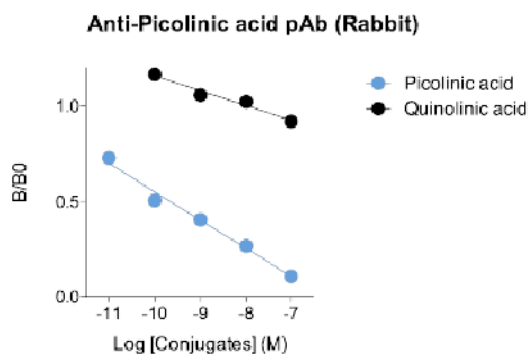
Product pictures

Picolinic acid



Picolinic acid

Affinity & specificity of Picolinic acid polyclonal antibody



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



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.