

Anti-conjugated L-Glutamate antibody

Ref: IS1001

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 L-Glutamate rabbit polyclonal antibody
Synonyms	Anti-L-Glutamic acid rabbit polyclonal antibody
Immunogen	Conjugated L-Glutamate
Specificity	When tested in competitive ELISA, the anti-conjugated L-Glutamate antibody did not show any significant cross reactivity with L-Glutamic acid analogs, including D-Glutamate, L-Glutamine and L-Aspartate 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

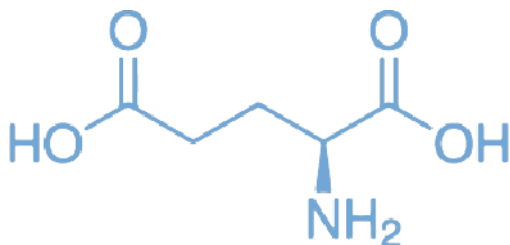
Antibody not yet cited.

Selected articles about L-Glutamate:

- Okubo Y et al. Imaging extrasynaptic glutamate dynamics in the brain. *Proc Natl Acad Sci U S A.* 2010 Apr 6;107(14):6526-31.
- Meldrum BS. Glutamate as a neurotransmitter in the brain: review of physiology and pathology. *J Nutr.* 2000 Apr;130(4S Suppl):1007S-15S.
- Hynd MR, Scott HL, Dodd PR. Glutamate-mediated excitotoxicity and neurodegeneration in Alzheimer's disease. *Neurochem Int.* 2004 Oct;45(5):583-95.

Pictures

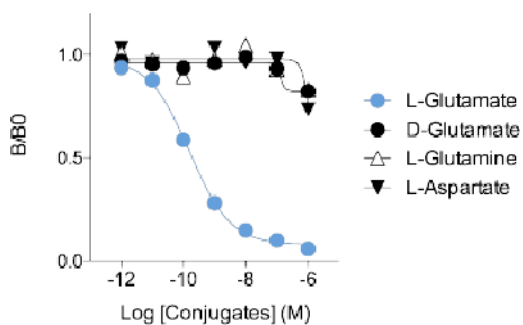
L-Glutamic acid (L-Glutamate)



L-Glutamate

Affinity & specificity of anti-L-Glutamate antibody

Anti-L-Glutamate pAb (Rabbit)



Competitive ELISA demonstrates that low amounts of L-Glutamate conjugate are required to abolish antigen-antibody reaction (high affinity), while rising concentrations of D-Glutamate, L-Glutamine and L-Aspartate 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.