

Anti-conjugated D-Glutamate pAb

Ref: IS1002

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

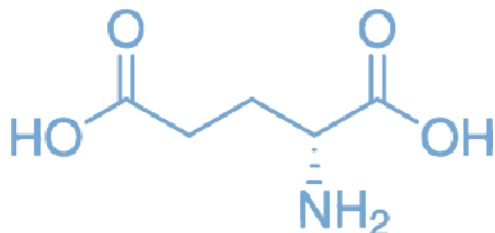
Antibody not yet cited.

Selected papers on D-Glutamic acid:

- [Dayer MR and Dayer MS. Hypoglycemic Effects of D-glutamic Acid in Diabetic Rats International Journal of Pharmacology, 2010 Nov 6: 930-936.](#)
- [Mangas A et al. Immunocytochemical visualization of D-glutamate in the rat brain. Neuroscience. 2007 Jan 19;144\(2\):654-64. Epub 2006 Nov 3.](#)
- [Fisher GH et al. Presence of D-aspartate and D-glutamate in tumor proteins. Cancer Biochem Biophys. 1995 Nov;15\(2\):79-82.](#)
- [Kera Y et al. Presence of free D-glutamate and D-aspartate in rat tissues. Biochim Biophys Acta. 1995 Feb 23;1243\(2\):283-6.](#)

Pictures

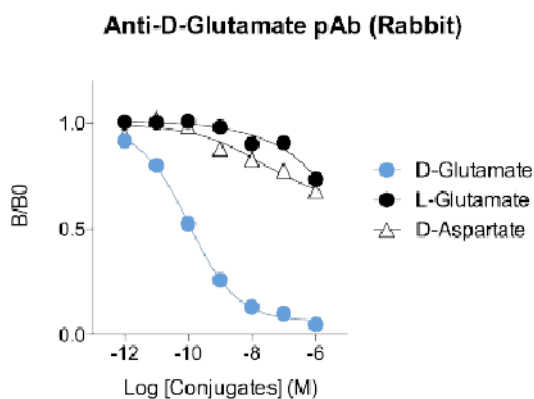
D-Glutamate (D-Glutamic acid)



D-Glutamate

Affinity & specificity of anti-D-Glutamate antibody

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



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.