

L-Kynurenine polyclonal antibody

Ref: IS1012

Validated for IHC in human tumor tissues, the anti-L-Kynurenine (Kyn) rabbit polyclonal antibody proved to work at 1/500 dilution on paraffin-embedded sections, a single vial thus catering for approximately 100 stainings.

| | |
|-----------------------------|--|
| Clonality | Polyclonal |
| Host | Rabbit (see anti-KYN mouse mAb) |
| Valided applications | <u>IHC</u> |
| Specie reactivity | Reacts with all species |
| References | Not yet cited to our knowledge. Submit content and <u>get a 10% discount!</u> |
| Format | 50µl |

Product information

Product overview

| | |
|---------------------|--|
| Product name | L-Kynurenine polyclonal antibody |
| Synonyms | KYN polyclonal antibody (S)-Kynurenine polyclonal antibody L-2-Amino-4-(2-aminophenyl)-4-oxobutanoic acid antibody Kynurenine polyclonal antibody ?-Anthraniloyl-L-alanine polyclonal antibody |
| Immunogen | Conjugated L-Kynurenine |
| Specificity | When tested in competitive ELISA, the anti-L-Kynurenine polyclonal antibody did not show any significant cross reactivity with competitors including anthranilic acid, 3-OH-Kynurenine conjugate and Tryptophan. |
| Lot number | 140401 |

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. |

Protocols

| | |
|-----------------------------------|--|
| Immunohistochemistry (IHC) | Dilute at 1:100-1:1000. Perform heat antigen retrieval (pH=6) before initiating IHC staining protocol on paraffin-embedded and frozen sections |
| Comments | Optimal working dilutions must be determined by the end-user |
| Restrictions | For research use only |

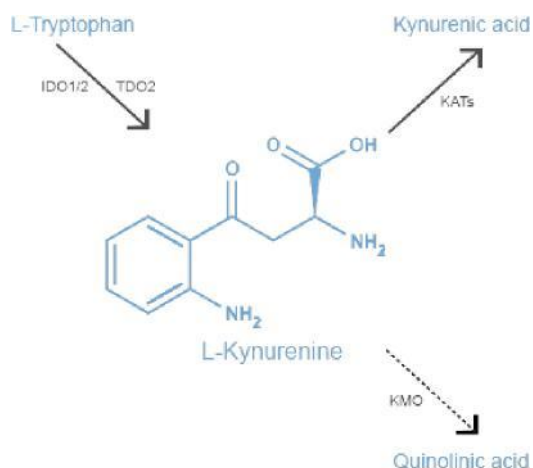
References

Antibody not yet cited. Submit an article and [get a 10% discount](#).

Selected articles on L-Kynurenine:

- [Munn DH, Mellor AL. Indoleamine 2,3 dioxygenase and metabolic control of immune responses. Trends Immunol. 2013 Mar;34\(3\):137-43. doi: 10.1016/j.it.2012.10.001. Epub 2012 Oct 25.](#)
- [Adams S, Braidly N, Bessede A, Brew BJ, Grant R, Teo C, Guillemin GJ. The kynurenine pathway in brain tumor pathogenesis. Cancer Res. 2012 Nov 15;72\(22\):5649-57. doi: 10.1158/0008-5472.CAN-12-0549. Epub 2012 Nov 9.](#)
- [Schwarcz R, Bruno JP, Muchowski PJ, Wu HQ. Kynurenines in the mammalian brain: when physiology meets pathology. Nat Rev Neurosci. 2012 Jul;13\(7\):465-77. doi: 10.1038/nrn3257.](#)
- [Opitz CA et al. An endogenous tumour-promoting ligand of the human aryl hydrocarbon receptor. Nature. 2011 Oct 5;478\(7368\):197-203. doi: 10.1038/nature10491.](#)
- [Favre D et al. Tryptophan catabolism by indoleamine 2,3-dioxygenase 1 alters the balance of TH17 to regulatory T cells in HIV disease. Sci Transl Med. 2010 May 19;2\(32\):32ra36. doi: 10.1126/scitranslmed.3000632.](#)

Product pictures

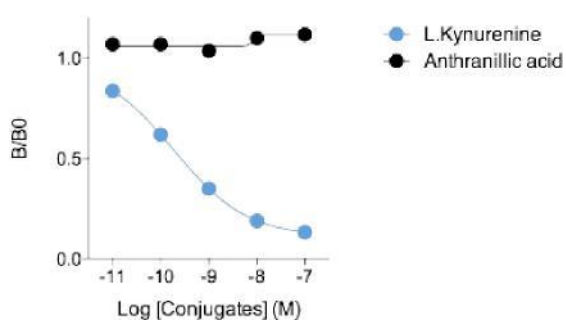


L-Kynurenine (KYN)

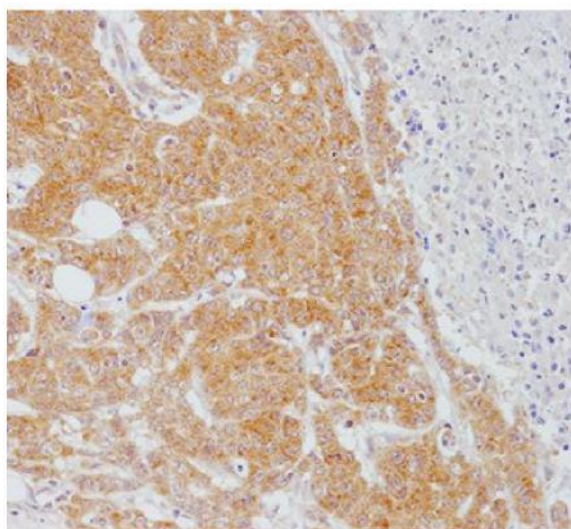
L-Kynurenine, the first stable metabolite of the kynurenine pathway, is synthesized from L-Tryptophan by indoleamine 2,3-dioxygenase (IDO1/2) or tryptophan 2,3-dioxygenase (TDO2) enzymes. Used as a biomarker of tryptophan degradation through the kynurenine pathway in various immune-related disorders, L-Kynurenine is known to activate the Aryl hydrocarbon Receptor (AhR), thereby exerting immunomodulatory effects and potentially promoting tumor immune escape. As a precursor of both neurotoxic and neuroactive metabolites, L-kynurenine is also monitored in the context of CNS disorders such as Parkinson's and Alzheimer's diseases, depression, schizophrenia, ...

Affinity & specificity of L-Kynurenine polyclonal antibody

Anti-L-Kynurenine pAb (Rabbit)



Competitive ELISA shows that low amounts of L-Kynurenine conjugate are required to abolish antigen-antibody reaction (high affinity), while rising concentrations of anthranilic acid conjugate do not affect reaction (high specificity).



L-Kynurenine accumulation in human colorectal tumor by IHC

Immunohistochemical analysis reveals cytoplasmic accumulation of L-Kynurenine in tumour cells in human colorectal cancer tissue. Paraffin-embedded tissue section was subjected to pH=6 antigen retrieval, and overnight incubation with primary anti-KYN polyclonal antibody (1/500 dilution). A polymer-conjugated secondary Ab was added and immunostaining was revealed using DAB.

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](mailto:info@eaglebio.com) or at 866-411-8023.