

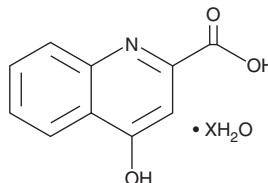
PRODUCT INFORMATION



Kynurenic Acid (hydrate)

Item No. 16792

CAS Registry No.: 345909-35-5
Formal Name: 4-hydroxy-2-quinolinecarboxylic acid, hydrate
Synonym: KYNA
MF: C₁₀H₇NO₃ • XH₂O
FW: 189.2
Purity: ≥98%
UV/Vis.: λ_{max}: 217, 247, 341 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Kynurenic acid (hydrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the kynurenic acid (hydrate) in the solvent of choice, which should be purged with an inert gas. Kynurenic acid (hydrate) is soluble in 0.1 M NaOH at a concentration of approximately 1 mg/ml.

Description

Kynurenic acid is an active metabolite of tryptophan (Item No. 29600 | 31210).¹ It is formed from tryptophan via a kynurenine (Item No. 11305) intermediate by kynurenine aminotransferases (KATs). Kynurenic acid is an antagonist of the NMDA and AMPA receptors as well as α7 nicotinic acetylcholine receptors (nAChRs; EC₅₀s = 235, 101, and 7 μM, respectively).² It is also an agonist of the aryl hydrocarbon receptor (AhR) and G protein-coupled receptor 35 (GPR35; EC₅₀s = 1.4 and 39 μM, respectively). It prevents lesioned hemisphere weight loss in a neonatal rat model of cerebral hypoxic-ischemia induced by carotid artery ligation when administered at a dose of 300 mg/kg.³ Kynurenic acid (1 and 5 mg/ml) is protective against rhabdomyer neurodegeneration in the eye in an *Htt*^{93Q} transgenic *Drosophila* model of Huntington's disease.⁴ Levels of kynurenic acid in cerebrospinal fluid are increased in patients with schizophrenia.⁵

References

- Schwarz, R., Bruno, J.P., Muchowski, P.J., *et al.* Kynurenines in the mammalian brain: When physiology meets pathology. *Nat. Rev. Neurosci.* **13**(7), 465-477 (2012).
- Turski, M.P., Turska, M., Paluszkiwicz, P., *et al.* Kynurenic acid in the digestive system-new facts, new challenges. *Int. J. Tryptophan Res.* **6**, 47-55 (2013).
- Andiné, P., Lehmann, A., Ellrén, K., *et al.* The excitatory amino acid antagonist kynurenic acid administered after hypoxic-ischemia in neonatal rats offers neuroprotection. *Neurosci. Lett.* **90**(1-2), 208-212 (1988).
- Campana, S., Green, E.W., Breda, C., *et al.* The kynurenine pathway modulates neurodegeneration in a *Drosophila* model of Huntington's disease. *Curr. Biol.* **21**(11), 9611-9666 (2011).
- Nilsson, L.K., Linderholm, K.R., Engberg, G., *et al.* Elevated levels of kynurenic acid in the cerebrospinal fluid of male patients with schizophrenia. *Schizophr. Res.* **80**(2-3), 315-322 (2005).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/02/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM