

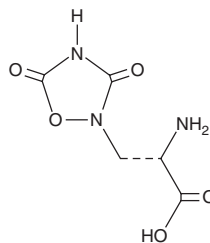
PRODUCT INFORMATION



L-Quisqualic Acid

Item No. 20211

CAS Registry No.: 52809-07-1
Formal Name: αS-amino-3,5-dioxo-1,2,4-oxadiazolidine-2-propanoic acid
MF: C₅H₇N₃O₅
FW: 189.1
Purity: ≥99%
Supplied as: A 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

L-Quisqualic acid is supplied as a solid. The solubility of L-quisqualic acid in water is approximately 25mM. We do not recommend storing the aqueous solution for more than one day.

Description

L-Quisqualic acid is a natural analog of glutamate that acts as an agonist at AMPA-selective and metabotropic glutamate receptors (EC₅₀s = 170, 10, 40, and 29 nM at GluR-A, mGluR1, mGluR3, and mGluR5, respectively), as well as the ionotropic kainate receptor (GRIK4; K_i = 6.43 nM).¹⁻⁶ L-Quisqualic acid is used to study receptor dynamics and as an excitotoxin to selectively destroy neurons in the brain or spinal cord.^{7,8}

References

1. Schoepp, D.D., Jane, D.E., and Monn, J.A. Pharmacological agents acting at subtypes of metabotropic glutamate receptors. *Neuropharmacology* **38(10)**, 1431-1476 (1999).
2. Bräuner-Osborne, H., Egebjerg, J., Nielsen, E.O., *et al.* Ligands for glutamate receptors: Design and therapeutic prospects. *J. Med. Chem.* **43(14)**, 2609-2645 (2000).
3. Bertrand, H.-O., Besis, A.-S., Pin, J.-P., *et al.* Common and selective molecular determinants involved in metabotropic glutamate receptor agonist activity. *J. Med. Chem.* **45(15)**, 3171-3183 (2002).
4. Kamboj, R.K., Schoepp, D.D., Nutt, S., *et al.* Molecular cloning, expression, and pharmacological characterization of humEAA1, a human kainate receptor subunit. *J. Neurochem.* **62(1)**, 1-9 (1994).
5. Mutel, V., Ellis, G.J., Adam, G., *et al.* Characterization of [³H]Quisqualate binding to recombinant rat metabotropic glutamate 1a and 5a receptors and to rat and human brain sections. *J. Neurochem.* **75(6)**, 2590-2601 (2000).
6. Uchino, S., Sakimura, K., Nagahari, K., *et al.* Mutations in a putative agonist binding region of the AMPA-selective glutamate receptor channel. *FEBS Lett.* **308(3)**, 253-257 (1992).
7. Lee, J.-W., Furmanski, O., Castellanos, D.A., *et al.* Prolonged nociceptive responses to hind paw formalin injection in rats with a spinal cord injury. *Neurosci. Lett.* **439(2)**, 212-215 (2008).
8. Muir, J.L., Page, K.J., Sirinathsinghji, D.J.S., *et al.* Excitotoxic lesions of basal forebrain cholinergic neurons: Effects on learning, memory and attention. *Behav. Brain Res.* **57(2)**, 123-131 (1993).

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, 12/19/2022

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