

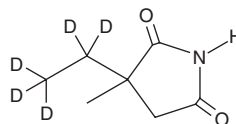
PRODUCT INFORMATION



Ethosuximide-d₅

Item No. 30759

CAS Registry No.: 1989660-59-4
Formal Name: 3-(ethyl-1,1,2,2,2-d₅)-3-methyl-2,5-pyrrolidinedione
Synonyms: CI-366-d₅, NSC 64013-d₅
MF: C₇H₆D₅NO₂
FW: 146.2
Chemical Purity: ≥98% (Ethosuximide)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
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

Ethosuximide-d₅ is intended for use as an internal standard for the quantification of ethosuximide (Item No. 23947) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Ethosuximide-d₅ is supplied as a solid. A stock solution may be made by dissolving the ethosuximide-d₅ in the solvent of choice, which should be purged with an inert gas. Ethosuximide-d₅ is soluble in methanol.

Description

Ethosuximide is an anticonvulsant.¹⁻⁵ It increases glucose, fructose-1,6-bisphosphate, and pyruvate levels in rat brain when administered at a dose of 200 mg/kg.¹ Ethosuximide (400 mg/kg) reduces the severity of audiogenic seizures in a rat model of barbiturate withdrawal-induced convulsions.² It also inhibits tonic hindlimb extension induced by pentylenetetrazole (PTZ; Item No. 18682) or brainstem stimulation (ED₅₀s = 35 and 70 mg/kg, respectively), as well as leptazol-induced clonic seizures (ED₅₀ = 230 mg/kg), in rats.^{3,4} Ethosuximide reduces resting tremor by 60% in a macaque model of Parkinson's disease induced by MPTP when administered at a dose of 150 mg/animal for 5 days.⁵ Formulations containing ethosuximide have been used in the treatment of petit mal seizures.

References

1. Nahorski, S.R. Biochemical effects of the anticonvulsants trimethadione, ethosuximide and chlorthalidoxepoxide in rat brain. *J. Neurochem.* **19(8)**, 1937-1946 (1972).
2. Norton, P.R. The effects of drugs on barbiturate withdrawal convulsions in the rat. *J. Pharm. Pharmacol.* **22(10)**, 763-766 (1970).
3. Consroe, P.F. and Wolkin, A.L. Anticonvulsant interaction of cannabidiol and ethosuximide in rats. *J. Pharm. Pharmacol.* **29(8)**, 500-501 (1977).
4. Chiu, P. and Burnham, W.M. The effect of anticonvulsant drugs on convulsions triggered by direct stimulation of the brainstem. *Neuropharmacology* **21(4)**, 355-359 (1982).
5. Gomez-Mancilla, B., Latulippe, J.F., Boucher, R., et al. Effect of ethosuximide on rest tremor in the MPTP monkey model. *Mov. Disord.* **7(2)**, 137-141 (1992).

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, 12/14/2022

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