

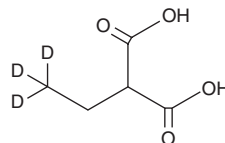
PRODUCT INFORMATION



Ethylmalonic Acid-d₃

Item No. 43923

CAS Registry No.: 70907-93-6
Formal Name: ethyl-2,2,2-d₃-propanedioic acid
Synonym: EMA-d₃
MF: C₅H₅D₃O₄
FW: 135.1
Chemical Purity: ≥98% (Ethylmalonic Acid)
Deuterium
Incorporation: ≥99% deuterated forms (d₁-d₃); ≤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

Ethylmalonic acid-d₃ is intended for use as an internal standard for the quantification of ethylmalonic acid 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).

Ethylmalonic acid-d₃ is supplied as a solid. A stock solution may be made by dissolving the ethylmalonic acid-d₃ in the solvent of choice, which should be purged with an inert gas. Ethylmalonic acid-d₃ is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

Ethylmalonic acid is a dicarboxylic acid and metabolite of the essential amino acid L-isoleucine and branched-chain amino acid L-alloisoleucine (Item No. 34904).¹ It is formed from L-isoleucine and L-alloisoleucine via an R-2-oxo-3-methylvaleric acid intermediate. Urinary levels of ethylmalonic acid are increased in patients with ethylmalonic encephalopathy, an inborn error of metabolism characterized by developmental delay, hypotonia, vascular instability, petechiae, acrocyanosis, chronic diarrhea, and lactic acidemia. Urinary levels of ethylmalonic acid are also increased in patients with short-chain acyl-CoA dehydrogenase deficiency.²

References

1. Nowaczyk, M.J.M., Lehotay, D.C., Platt, B.-A., *et al.* Ethylmalonic and methylsuccinic aciduria in ethylmalonic encephalopathy arise from abnormal isoleucine metabolism. *Metabolism* **47(7)**, 836-839 (1998).
2. Gallant, N.M., Leydiker, K., Tang, H., *et al.* Biochemical, molecular, and clinical characteristics of children with short chain acyl-CoA dehydrogenase deficiency detected by newborn screening in California. *Mol. Genet. Metab.* **106(1)**, 55-61 (2012).

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, 07/16/2025

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