

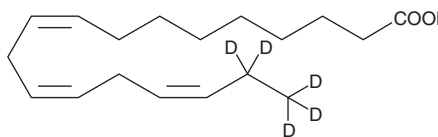
PRODUCT INFORMATION



α -Linolenic Acid-d₅

Item No. 9002286

CAS Registry No.: 145191-04-4
Formal Name: 9Z,12Z,15Z-octadecatrienoic-17,17',18,18,18-d₅ acid
Synonyms: ALA-d₅, FA 18:3-d₅
MF: C₁₈H₂₅D₅O₂
FW: 283.5
Chemical Purity: ≥98% α -Linolenic Acid
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

α -Linolenic acid-d₅ (ALA-d₅) is intended for use as an internal standard for the quantification of ALA (Item No. 90210) 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).

ALA-d₅ is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of ALA-d₅ in these solvents is approximately 100 mg/ml.

Description

ALA is an essential fatty acid found in leafy green vegetables. ALA, as part of a low saturated fat diet, helps prevent cardiovascular disease. ALA decreases blood pressure, serum cholesterol levels, and platelet aggregation.¹

Reference

1. Allman, M.A., Pena, M.M., Pang, D. Supplementation with flaxseed oil versus sunflowerseed oil in healthy young men consuming a low fat diet: Effects on platelet composition and function. *Eur. J. Clin. Nutr.* **49(3)**, 169-178 (1995).

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, 02/20/2024

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