

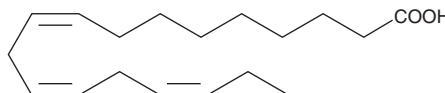
PRODUCT INFORMATION



α -Linolenic Acid

Item No. 90210

CAS Registry No.: 463-40-1
Formal Name: 9Z,12Z,15Z-octadecatrienoic acid
Synonyms: ALA, C18:3 n-3, C18:3 (9Z,12Z,15Z)
MF: C₁₈H₃₀O₂
FW: 278.4
Purity: \geq 98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

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

ALA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of ALA should be diluted with the aqueous buffer of choice. The solubility of ALA in PBS (pH 7.2) is approximately 100 μ g/ml and approximately 1 mg/ml in 0.15 M Tris-HCl (pH 8.5). We do not recommend storing the aqueous solution for more than one day. We do not recommend storing the aqueous solution for more than one day.

Description

ALA is an essential ω -3 polyunsaturated fatty acid found in plants.¹ It is converted to the longer-chain fatty acids eicosapentaenoic (Item Nos. 90110 | 90110.1 | 21908), docosapentaenoic (Item Nos. 90165 | 21907), and docosahexaenoic acid (Item No. 90310).² It is also catabolized *via* fatty acid β -oxidation for energy or the synthesis of saturated and monounsaturated fatty acids, or stored in adipose tissue. Dietary consumption of α -linolenic acid-containing foods is positively associated with a moderately lower risk of cardiovascular disease.¹

References

1. Pan, A., Chen, M., Chowdhury, R., *et al.* α -Linolenic acid and risk of cardiovascular disease: A systematic review and meta-analysis. *Am. J. Clin. Nutr.* **96**(6), 1262-1273 (2012).
2. Burdge, G.C. Metabolism of α -linolenic acid in humans. *Prostaglandins Leukot. Essent. Fatty Acids* **75**(3), 161-168 (2006).

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, 09/19/2023

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