

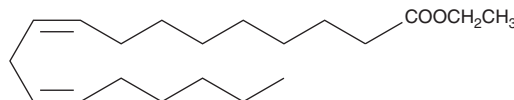
PRODUCT INFORMATION



Linoleic Acid ethyl ester

Item No. 10008198

CAS Registry No.: 544-35-4
Formal Name: 9Z,12Z-octadecadienoic acid, ethyl ester
Synonyms: C18:2 (cis-9,12) ethyl ester, Ethyl cis-9,12-Octadecadienoate, Ethyl Linoleate; Telfairic Acid ethyl ester, SFE 20:2
MF: C₂₀H₃₆O₂
FW: 308.5
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Linoleic acid ethyl ester 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 ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of linoleic acid ethyl ester in these solvents is approximately 100 mg/ml.

Description

Linoleic acid ethyl ester is an esterified form of linoleic acid (Item Nos. 90150 | 90150.1 | 21909). It is active against the bacteria *S. mutans*, *A. actinomycetemcomitans*, *P. gingivalis*, *S. gordonii*, and *S. sanguis* and the fungus *C. albicans* when used at a concentration of 25 µg/ml.¹ Linoleic acid ethyl ester (3 mg/kg) prevents arteriosclerosis in rabbits fed a high-cholesterol diet.²

References

1. Huang, C.B., George, B., and Ebersole, J.L. Antimicrobial activity of n-6, n-7 and n-9 fatty acids and their esters for oral microorganisms. *Arch. Oral Biol.* **55(8)**, 555-560 (2010).
2. Datey, K.K., Purandare, N.M., and Dalvi, C.P. Ethyl linoleate as an antiatherogenic agent. An experimental study in rabbits. *Angiology* **17(10)**, 732-737 (1966).

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, 03/27/2024

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