

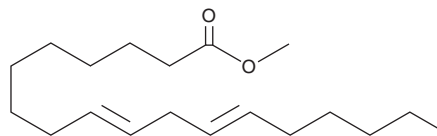
PRODUCT INFORMATION



Linoelaidic Acid methyl ester

Item No. 26732

CAS Registry No.: 2566-97-4
Formal Name: 9E,12E-octadecadienoic acid, methyl ester
Synonyms: C18:2 (*trans*-9,*trans*-12) methyl ester,
Methyl Linoelaidate,
Methyl *trans*-9,*trans*-12-Octadecadienoate,
SFE 19:2
MF: C₁₉H₃₄O₂
FW: 294.5
Purity: ≥95%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Linoelaidic acid methyl ester is supplied as a neat oil. A stock solution may be made by dissolving the linoelaidic acid methyl ester in the solvent of choice, which should be purged with an inert gas. Linoelaidic acid methyl ester is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of linoelaidic acid methyl ester in ethanol is approximately 50 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

Linoelaidic acid methyl ester is an esterified form of linoelaidic acid (Item No. 90160). It has been found in *H. sabdariffa* essential oil as well as biodiesel produced from sea mango seed oil.^{1,2}

References

1. Shen, C.Y., Zhang, T.T., Zhang, W.L., *et al.* Anti-inflammatory activities of essential oil isolated from the calyx of *Hibiscus sabdariffa* L. *Food Funct.* **7**(10), 4451-4459 (2016).
2. Lie, J., Rizkiana, B., Soetaredjo, F.E., *et al.* Production of biodiesel from sea mango (*Cerbera odollam*) seed using in situ subcritical methanol-water under a non catalytic process. *Int. J. Indust. Chem.* **9**(1), 53-59 (2018).

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, 03/04/2024

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