

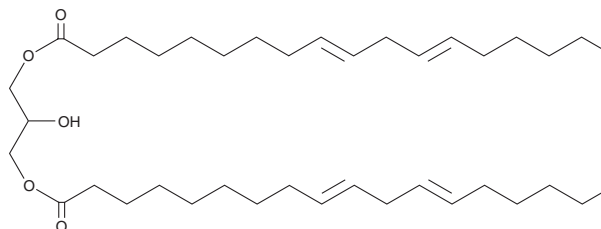
PRODUCT INFORMATION



1,3-Dilinoelaidoyl Glycerol

Item No. 26983

CAS Registry No.: 372490-73-8
Formal Name: (9E,9'E,12E,12'E)-9,12-octadecadienoic acid, 1,1'-(2-hydroxy-1,3-propanediyl) ester
Synonyms: DG(18:2/0:0/18:2), 1,3-Dilinoelaidin
MF: C₃₉H₆₈O₅
FW: 617.0
Purity: ≥98%
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

1,3-Dilinoelaidoyl glycerol is supplied as a solid. A stock solution may be made by dissolving the 1,3-dilinoelaidoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Dilinoelaidoyl glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 1,3-dilinoelaidoyl glycerol in these solvents is approximately 10 mg/ml.

1,3-Dilinoelaidoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,3-dilinoelaidoyl glycerol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 1,3-Dilinoelaidoyl glycerol has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

1,3-Dilinoelaidoyl glycerol is a diacylglycerol that contains linoelaidic acid (Item No. 90160) at the *sn*-1 and *sn*-3 positions. It is a collision-induced decomposition product of synthetic linoleic ammoniated triglycerides.¹

Reference

1. Gakwaya, R., Li, X., Wong, Y.L., *et al.* Examining the collision-induced decomposition spectra of ammoniated triglycerides. III. The linoleate and arachidonate series. *Rapid Commun. Mass Spectrom.* **21(20)**, 3262-3268 (2007).

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, 12/08/2022

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