

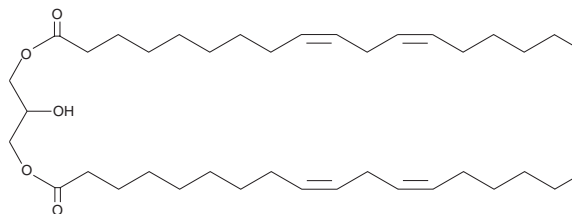
PRODUCT INFORMATION



1,3-Dilinoleoyl Glycerol

Item No. 16481

CAS Registry No.: 15818-46-9
Formal Name: 9Z,12Z-octadecadienoic acid, 1,1'-(2-hydroxy-1,3-propanediyl) ester
Synonyms: 1,3-Dilinolein, 1,3-Glycerol dilinoleate
MF: C₃₉H₆₈O₅
FW: 617.0
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

1,3-Dilinoleoyl glycerol is supplied as a solution in ethanol. To change the solvent, simply evaporate the 1,3-dilinoleoyl glycerol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol and dimethyl formamide purged with an inert gas can be used. The solubility of 1,3-dilinoleoyl glycerol in these solvents is approximately 10 mg/ml.

1,3-Dilinoleoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 1,3-dilinoleoyl glycerol should be diluted with the aqueous buffer of choice. 1,3-Dilinoleoyl 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-Dilinoleoyl glycerol is an unsaturated diacylglycerol with linoleic acid (Item No. 90150) (18:2) side-chains attached at both the C-1 and C-3 positions that is useful for lipid research.

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, 05/05/2023

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