

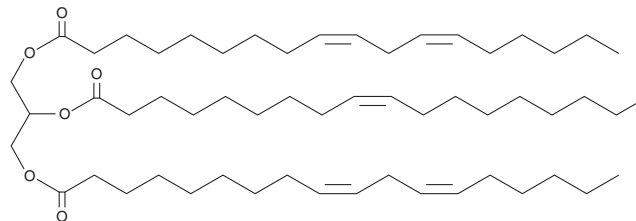
PRODUCT INFORMATION



1,3-Dilinoleoyl-2-Oleoyl Glycerol

Item No. 27064

CAS Registry No.: 2190-22-9
Formal Name: 9Z,12Z-octadecadienoic acid, 1,1'-[2-[[[(9Z)-1-oxo-9-octadecenyl]oxy]-1,3-propanediyl] ester
Synonyms: 1,3-Linolein-2-Olein, TG(18:2/18:1/18:2)
MF: C₅₇H₁₀₀O₆
FW: 881.4
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

1,3-Dilinoleoyl-2-oleoyl glycerol is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. 1,3-Dilinoleoyl-2-oleoyl glycerol is soluble in organic solvents such as ethanol and dimethyl formamide, which should be purged with an inert gas. The solubility of 1,3-dilinoleoyl-2-oleoyl glycerol in these solvents is approximately 10 mg/ml.

Description

1,3-Dilinoleoyl-2-oleoyl glycerol is a triacylglycerol that contains linoleic acid (Item Nos. 90150 | 90150.1 | 21909) at the *sn*-1 and *sn*-3 positions and oleic acid (Item Nos. 90260 | 24659) at the *sn*-2 position. It has been found in grape seed, pumpkin seed, soybean, sunflower, and wheat germ oils, with the highest content in grape seed oil.¹

Reference

1. Jakab, A., Jablonkai, I., and Forgács, E. Quantification of the ratio of positional isomer dilinoleoyl-oleoyl glycerols in vegetable oils. *Rapid Commun. Mass Spectrom.* **17(20)**, 2295-2302 (2003).

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, 12/07/2023

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