

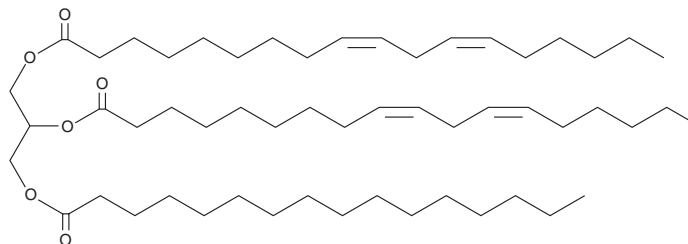
PRODUCT INFORMATION



1,2-Dilinoleoyl-3-Palmitoyl-*rac*-glycerol

Item No. 26887

CAS Registry No.: 2190-15-0
Formal Name: 9Z,12Z-octadecadienoic acid, 1,1'-[1-[[[1-oxohexadecyl)oxy] methyl]-1,2-ethanediyl] ester
Synonyms: 1,2-Linolein-3-Palmitin, TG(18:2/18:2/16:0)
MF: C₅₅H₉₈O₆
FW: 855.4
Purity: ≥95%
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

1,2-Dilinoleoyl-3-palmitoyl-*rac*-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. 1,2-Dilinoleoyl-3-palmitoyl-*rac*-glycerol is slightly soluble in chloroform and methanol.

Description

1,2-Dilinoleoyl-3-palmitoyl-*rac*-glycerol is a triacylglycerol that contains linoleic acid (Item Nos. 90150 | 90150.1 | 21909) at the *sn*-1 and *sn*-2 positions and palmitic acid (Item No. 10006627) at the *sn*-3 position. It has been found in a variety of vegetable oils, including poppy seed, hazelnut, maize, and olive oils.¹ 1,2-Dilinoleoyl-3-palmitoyl-*rac*-glycerol (3% w/v) reduces scald development on apples of the Delicious variety when applied immediately following harvest and assessed after six months of storage.²

References

- Mottram, H.R., Woodbury, S.E., and Evershed, R.P. Identification of triacylglycerol positional isomers present in vegetable oils by high performance liquid chromatography/atmospheric pressure chemical ionization mass spectrometry. *Rapid Commun. Mass Spectrom.* **11(12)**, 1240-1252 (1998).
- Ju, Z., Duan, Y., and Ju, Z. Mono-, di-, and tri-acylglycerols and phospholipids from plant oils inhibit scald development in 'Delicious' apples. *Postharvest Bio. Technol.* **19(1)** (2000).

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, 10/24/2022

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