

PRODUCT INFORMATION

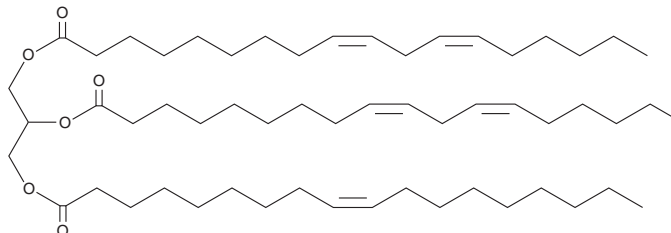


1,2-Dilinoleoyl-3-Oleoyl-*rac*-glycerol

Item No. 28178

CAS Registry No.: 2190-21-8
Formal Name: (9Z,12Z)-9,12-octadecadienoic acid, 1,1'-[1-[[[(9Z)-1-oxo-9-octadecen-1-yl]oxy]methyl]-1,2-ethanediyl] ester
Synonyms: 1,2-Dilinolein-3-Olein, TG(18:2/18:2/18:1)

MF: C₅₇H₁₀₀O₆
FW: 881.4
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-oleoyl-*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. A stock solution may be made by dissolving the 1,2-dilinoleoyl-3-oleoyl-*rac*-glycerol in the solvent of choice. 1,2-Dilinoleoyl-3-oleoyl-*rac*-glycerol is slightly soluble in methanol and water.

Description

1,2-Dilinoleoyl-3-oleoyl-*rac*-glycerol is a triacylglycerol that contains linoleic acid (Item Nos. 90150 | 90150.1 | 21909) at the *sn*-1 and *sn*-2 positions and oleic acid (Item Nos. 90260 | 24659) at the *sn*-3 position. It has been found in ostrich and emu oils, as well as in the fat body of male *B. lapidarius* bumblebees.^{1,2} 1,2-Dilinoleoyl-3-oleoyl-*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.³

References

1. Zhou, Y., Xue, Y., Chen, G.C., *et al.* Rapid separation and characterisation of triacylglycerols in ostrich oil by ultra performance liquid chromatography coupled with quadrupole time-of-flight mass spectrometry. *Food Chem.* **141**(3), 2098-2102 (2013).
2. Cvačka, J., Hovorka, O., Jiroš, P., *et al.* Analysis of triacylglycerols in fat body of bumblebees by chromatographic methods. *J. Chromatogr. A.* **1101**(1-2), 226-237 (2006).
3. 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), 1-7 (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, 03/22/2022

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