

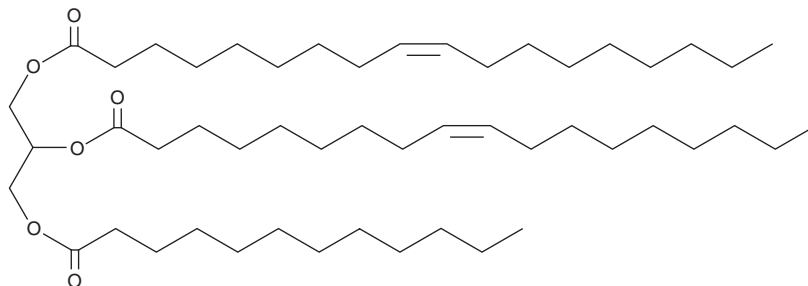
PRODUCT INFORMATION



1,2-Dioleoyl-3-Lauroyl-*rac*-glycerol

Item No. 28548

CAS Registry No.: 4016-52-8
Formal Name: 9Z-octadecenoic acid
1,1'-[1-[[[(1-oxododecyl)oxy]
methyl]-1,2-ethanediyl] ester
Synonyms: 1,2-Olein-3-Laurin,
TG(18:1/18:1/12:0)
MF: C₅₁H₉₄O₆
FW: 803.3
Purity: ≥98%
Supplied as: An oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

1,2-Dioleoyl-3-lauroyl-*rac*-glycerol is supplied as an oil. A stock solution may be made by dissolving the 1,2-dioleoyl-3-lauroyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1,2-Dioleoyl-3-lauroyl-*rac*-glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 1,2-dioleoyl-3-lauroyl-*rac*-glycerol in these solvents is approximately 10 mg/ml. 1,2-Dioleoyl-3-lauroyl-*rac*-glycerol is slightly soluble in chloroform.

1,2-Dioleoyl-3-lauroyl-*rac*-glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,2-dioleoyl-3-lauroyl-*rac*-glycerol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 1,2-Dioleoyl-3-lauroyl-*rac*-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

1,2-Dioleoyl-3-lauroyl-*rac*-glycerol is a triacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 and *sn*-2 positions and lauric acid (Item No. 10006626) at the *sn*-3 position. It has been found in date seed oil and the fat body of male *B. lapidarius* bumblebees.^{1,2}

References

- Holcapek, M., Lísá, M., Jandera, P., *et al.* Quantitation of triacylglycerols in plant oils using HPLC with APCI-MS, evaporative light-scattering, and UV detection. *J. Sep. Sci.* **28(12)**, 1315-1333 (2005).
- 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).

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, 10/27/2022

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