

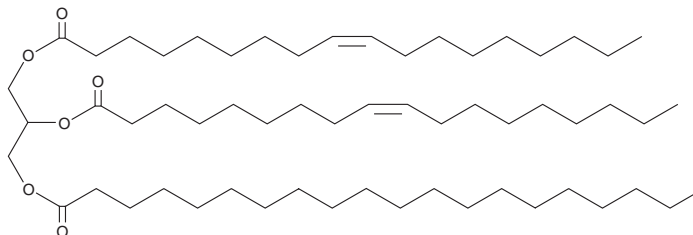
PRODUCT INFORMATION



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

Item No. 29060

CAS Registry No.: 77145-65-4
Formal Name: eicosanoic acid, 2,3-bis[[[9Z]-1-oxo-9-octadecen-1-yl]oxy]propyl ester
Synonyms: 1,2-Olein-3-Arachidin, TG(18:1/18:1/20:0)
MF: C₅₉H₁₁₀O₆
FW: 915.5
Purity: ≥95%
Supplied as: A crystalline solid
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-arachidoyl-*rac*-glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-dioleoyl-3-arachidoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1,2-Dioleoyl-3-arachidoyl-*rac*-glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 1,2-dioleoyl-3-arachidoyl-*rac*-glycerol in these solvents is approximately 10 mg/ml. It is also slightly soluble in chloroform.

1,2-Dioleoyl-3-arachidoyl-*rac*-glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,2-dioleoyl-3-arachidoyl-*rac*-glycerol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 1,2-Dioleoyl-3-arachidoyl-*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-arachidoyl-*rac*-glycerol is a triacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 and *sn*-2 positions and arachidic acid (Item Nos. 9000339 | 21906) at the *sn*-3 position. It has been found in hazelnut, cashew, date seed, corn, sunflower, and soybean oils.^{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).
- Gao, B., Luo, Y., Lu, W., *et al.* Triacylglycerol compositions of sunflower, corn and soybean oils examined with supercritical CO₂ ultra-performance convergence chromatography combined with quadrupole time-of-flight mass spectrometry. *Food Chem.* **218**, 569-574 (2017).

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, 10/28/2022

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