

PRODUCT INFORMATION



1-Arachidoyl-2-Oleoyl-3-Palmitoyl-*rac*-glycerol

Item No. 28556

CAS Registry No.: 81655-42-7
Formal Name: eicosanoic acid
3-[[[1-oxohexadecyl]oxy]-2-[[[(9Z)-1-oxo-9-octadecen-1-yl]oxy]propyl ester

Synonyms: 1-Arachidin-2-Olein-3-Palmitin, TG(20:0/18:1/16:0)

MF: C₅₇H₁₀₈O₆

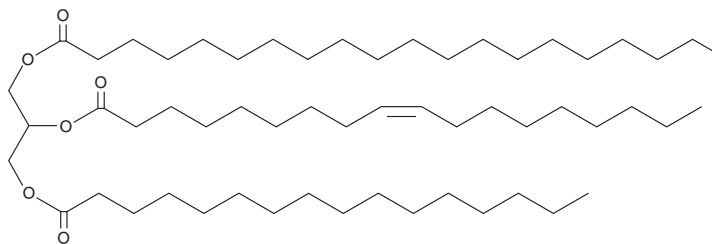
FW: 889.5

Purity: ≥98%

Supplied as: A 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-Arachidoyl-2-oleoyl-3-palmitoyl-*rac*-glycerol is supplied as a solid. A stock solution may be made by dissolving the 1-arachidoyl-2-oleoyl-3-palmitoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1-Arachidoyl-2-oleoyl-3-palmitoyl-*rac*-glycerol is soluble in chloroform and methanol.

Description

1-Arachidoyl-2-oleoyl-3-palmitoyl-*rac*-glycerol is a triacylglycerol that contains arachidic acid (Item Nos. 9000339 | 21906), oleic acid (Item Nos. 90260 | 24659), and palmitic acid (Item No. 10006627) at the *sn*-1, *sn*-2, and *sn*-3 positions, respectively. It has been found in a variety of plant oils, including poppy seed, hazelnut, maize, olive, yellow melon, and date seed oils.^{1,2}

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 (1997).
- Holcapek, M., Lisa, 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).

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, 12/05/2022

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