

PRODUCT INFORMATION



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

Item No. 28553

CAS Registry No.: 669762-61-2
Formal Name: eicosanoic acid, 1-[[[(1-oxohexadecyl)oxy]methyl]-2-[[[(9Z)-1-oxo-9-octadecen-1-yl]oxy]ethyl] ester

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

MF: C₅₇H₁₀₈O₆

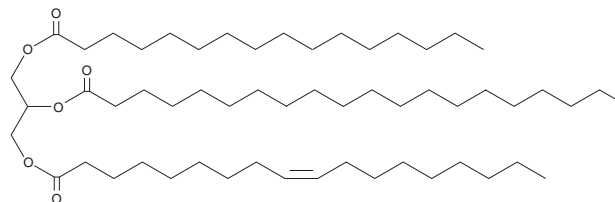
FW: 889.5

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-Palmitoyl-2-arachidoyl-3-oleoyl-*rac*-glycerol is supplied as a solid. A stock solution may be made by dissolving the 1-palmitoyl-2-arachidoyl-3-oleoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1-Palmitoyl-2-arachidoyl-3-oleoyl-*rac*-glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 1-palmitoyl-2-arachidoyl-3-oleoyl-*rac*-glycerol in these solvents is approximately 2 mg/ml. 1-Palmitoyl-2-arachidoyl-3-oleoyl-*rac*-glycerol is also slightly soluble in DMSO.

Description

1-Palmitoyl-2-arachidoyl-3-oleoyl-*rac*-glycerol is a triacylglycerol that contains palmitic acid (Item No. 10006627), arachidic acid (Item No. 9000339 | 21906), and oleic acid (Item Nos. 90260 | 24659) at the *sn*-1, *sn*-2, and *sn*-3 positions, respectively. It has been found in cod liver oil.¹

Reference

1. Zeng, Y.X., Araujo, P., Du, Z.Y., *et al.* Elucidation of triacylglycerols in cod liver oil by liquid chromatography electrospray tandem ion-trap mass spectrometry. *Talanta* **82(4)**, 1261-1270 (2010).

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, 11/04/2022

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