

PRODUCT INFORMATION



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

Item No. 29057

CAS Registry No.: 675876-04-7
Formal Name: 5Z,8Z,11Z,14Z-eicosatetraenoic acid, 3-[(1-oxohexadecyl)oxy]-2-[[[(9Z)-1-oxo-9-octadecen-1-yl]oxy]propyl ester

Synonyms: 1-Palmitin-2-Olein-3-Arachidonin, TG(16:0/18:1/20:4)

MF: C₅₇H₁₀₀O₆

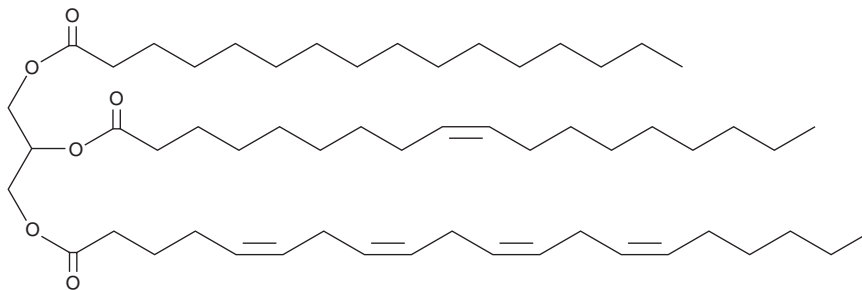
FW: 881.4

Purity: ≥95%

Supplied as: A solution in ethanol

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-oleoyl-3-arachidonoyl-*rac*-glycerol is supplied as a solution in ethanol. To change the solvent, simply evaporate the XX under a gentle stream of nitrogen and immediately add the solvent of choice. 1-Palmitoyl-2-oleoyl-3-arachidonoyl-*rac*-glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 1-palmitoyl-2-oleoyl-3-arachidonoyl-*rac*-glycerol in these solvents is approximately 10 mg/ml. 1-Palmitoyl-2-oleoyl-3-arachidonoyl-*rac*-glycerol is slightly soluble in chloroform.

1-Palmitoyl-2-oleoyl-3-arachidonoyl-*rac*-glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1-palmitoyl-2-oleoyl-3-arachidonoyl-*rac*-glycerol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 1-Palmitoyl-2-oleoyl-3-arachidonoyl-*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-Palmitoyl-2-oleoyl-3-arachidonoyl-*rac*-glycerol is a triacylglycerol that contains palmitic acid (Item No. 10006627), oleic acid (Item Nos. 90260 | 24659), and arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) at the *sn*-1, *sn*-2, and *sn*-3 positions, respectively. It has been found in lipid extracts from whole rat retina.¹

Reference

1. Armstrong, D. *Lipidomics Volume 1: Methods and Protocols*. 1st ed., Humana Press, New York, NY, (2009).

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, 02/20/2024

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