

PRODUCT INFORMATION



1,3-Dipalmitoyl-2-Arachidonoyl-*sn*-glycerol

Item No. 29341

Formal Name: 2-(((5Z,8Z,11Z,14Z)-icosa-5,8,11,14-tetraenyl)oxy)propane-

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

MF: C₅₅H₉₈O₆

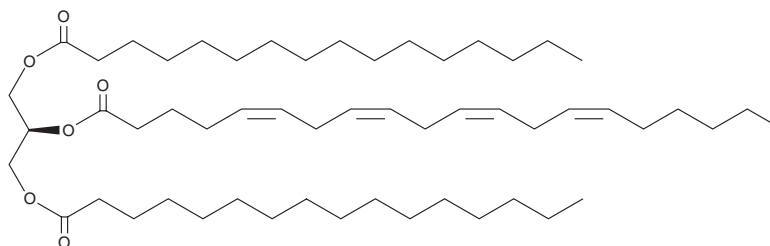
FW: 855.4

Purity: ≥98%

Supplied as: A solution in methyl acetate

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

1,3-Dipalmitoyl-2-arachidonoyl-*sn*-glycerol is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of 1,3-dipalmitoyl-2-arachidonoyl-*sn*-glycerol in DMF is approximately 10 mg/ml.

Description

1,3-Dipalmitoyl-2-arachidonoyl-*sn*-glycerol is a triacylglycerol that contains palmitic acid (Item No. 10006627) at the *sn*-1 and *sn*-3 positions and arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) at the *sn*-2 position.

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, 11/09/2023

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