

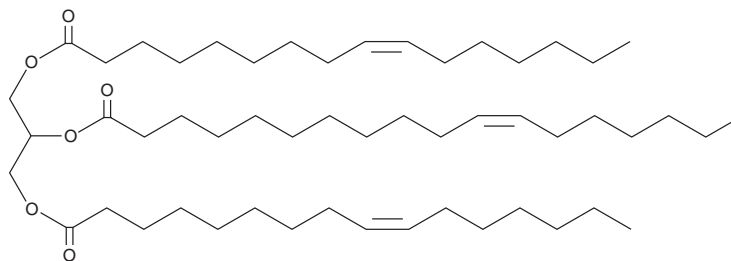
PRODUCT INFORMATION



1,3-Dipalmitoleoyl-2-11(Z)-Octadecenoyl Glycerol

Item No. 27380

CAS Registry No.: 2348469-98-5
Formal Name: (11Z)-11-octadecenoic acid, 2-[[[(9Z)-1-oxo-9-hexadecen-1-yl]oxy]-1-[[[(9Z)-1-oxo-9-hexadecen-1-yl]oxy]methyl] ethyl ester
Synonyms: 1,3-Palmitolein-2-Vaccenate, TG(16:1/18:1/16:1)
MF: C₅₃H₉₆O₆
FW: 829.3
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

1,3-Dipalmitoleoyl-2-11(Z)-octadecenoyl 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 ethanol and dimethyl formamide purged with an inert gas can be used. The solubility of 1,3-dipalmitoleoyl-2-11(Z)-octadecenoyl glycerol in these solvents is approximately 10 mg/ml. 1,3-Dipalmitoleoyl-2-11(Z)-octadecenoyl glycerol is slightly soluble in chloroform.

1,3-Dipalmitoleoyl-2-11(Z)-octadecenoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the solution of 1,3-dipalmitoleoyl-2-11(Z)-octadecenoyl glycerol should be diluted with the aqueous buffer of choice. 1,3-Dipalmitoleoyl-2-11(Z)-octadecenoyl glycerol has a solubility of 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

1,3-Dipalmitoleoyl-2-11(Z)-octadecenoyl glycerol is a triacylglycerol that contains palmitoleic acid (Item Nos. 10009871 | 21911) at the *sn*-1 and *sn*-3 positions and *cis*-vaccenic acid (Item No. 20023) 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, 12/14/2023

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