

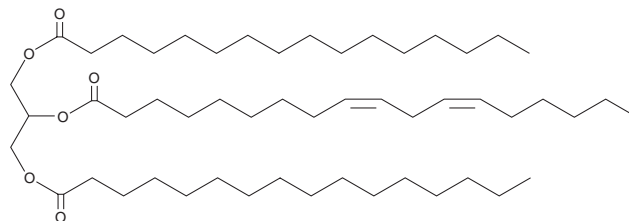
PRODUCT INFORMATION



1,3-Dipalmitoyl-2-Linoleoyl Glycerol

Item No. 27072

CAS Registry No.: 2442-56-0
Formal Name: 9Z,12Z-octadecadienoic acid, 2-[[1-oxohexadecyl)oxy]-1-[[1-oxohexadecyl)oxy]methyl]ethyl ester
Synonyms: 1,3-Palmitin-2-Linolein, TG(16:0/18:2/16:0)
MF: C₅₃H₉₈O₆
FW: 831.3
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-linoleoyl 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. 1,3-Dipalmitoyl-2-linoleoyl glycerol is slightly soluble in chloroform and methanol.

Description

1,3-Dipalmitoyl-2-linoleoyl glycerol is a triacylglycerol that contains palmitic acid (Item No. 10006627) at the *sn*-1 and *sn*-3 positions and linoleic acid (Item Nos. 90150 | 90150.1 | 21909) at the *sn*-2 position. It has been found in the palm stearin fraction of palm oil and as a minor component of soybean germ oil.^{1,2}

References

1. Dian, N.L.H.M., Lin, S.W., Let, C.C., *et al.* Characteristics of simple and chemically interesterified blends containing palm stearin, sunflower oil and palm kernel olein and potential application of the blends in fats spread formulation. *Oil Palm Bulletin* **43**, 38-49 (2018).
2. Yu, D., Elfalleh, W., He, S., *et al.* Physicochemical properties and minor lipid components of soybean germ. *J. Am. Oil. Chem. Soc.* **90(10)**, 1551-1558 (2013).

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, 10/13/2022

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