

PRODUCT INFORMATION



1,3-Dioleoyl-2-Linoleoyl Glycerol

Item No. 26972

CAS Registry No.: 2190-19-4
Formal Name: 9Z,12Z-octadecadienoic acid, 2-[[[(9Z)-1-oxo-9-octadecen-1-yl]oxy]-1-[[[(9Z)-1-oxo-9-octadecen-1-yl]oxy]methyl]ethyl ester

Synonyms: 1,3-Olein-2-Linolein, OLO, TG(18:1/18:2/18:1), 18:1/18:2/18:1-TG

MF: C₅₇H₁₀₂O₆

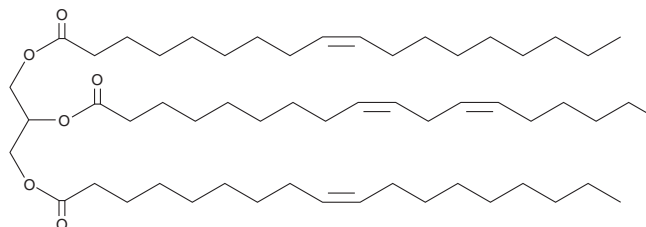
FW: 883.4

Purity: ≥95%

Supplied as: A solution in hexane

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-Dioleoyl-2-linoleoyl glycerol is supplied as a solution in hexane. To change the solvent, simply evaporate the hexane under a gentle stream of nitrogen and immediately add the solvent of choice. 1,3-Dioleoyl-2-linoleoyl glycerol is sparingly soluble (1-10 mg/ml) in chloroform.

Description

1,3-Dioleoyl-2-linoleoyl glycerol is a triacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) 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 as a major molecular species in butterfat.¹ It has also been found in *A. trifoliata* seed oil (ASO), as well as, to a lesser extent, in physical blends and interesterified products of ASO.²

References

1. Kalo, P., Kemppinen, A., and Ollilainen, V. Determination of triacylglycerols in butterfat by normal-phase HPLC and electrospray-tandem mass spectrometry. *Lipids* **44**(2), 169-195 (2009).
2. Zhao, S.-Q., Hu, J.-N., Zhu, X.-M., *et al.* Characteristics and feasibility of trans-free plastic fats through lipozyme TL IM-catalyzed interesterification of palm stearin and *Akebia trifoliata* variety australis seed oil. *J. Agric. Food Chem.* **62**(14), 3293-3300 (2014).

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, 03/18/2024

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