

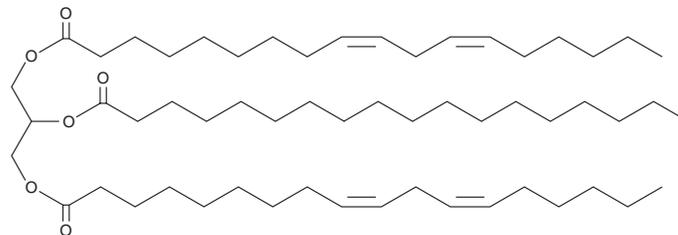
PRODUCT INFORMATION



1,3-Dilinoleoyl-2-Stearoyl Glycerol

Item No. 26971

CAS Registry No.: 2190-18-3
Formal Name: 9,12-octadecadienoic acid, 1,1'-[2-[(1-oxooctadecyl)oxy]-1,3-propanediyl] ester
Synonyms: 1,3-Linolein-2-Stearin, 18:2/18:0/18:2-TG, TG(18:2/18:0/18:2)
MF: C₅₇H₁₀₂O₆
FW: 883.4
Purity: ≥98%
Supplied as: A solution in ethanol
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-Dilinoleoyl-2-stearoyl glycerol is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice, which should be purged with an inert gas. 1,3-Dilinoleoyl-2-stearoyl glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 1,3-dilinoleoyl-2-stearoyl glycerol in these solvents is approximately 10 mg/ml. 1,3-Dilinoleoyl-2-stearoyl glycerol is also slightly soluble in chloroform.

1,3-Dilinoleoyl-2-stearoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 1,3-dilinoleoyl-2-stearoyl glycerol should be diluted with the aqueous buffer of choice. 1,3-Dilinoleoyl-2-stearoyl 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-Dilinoleoyl-2-stearoyl glycerol is a triacylglycerol that contains linoleic acid (Item Nos. 90150 | 90150.1 | 21909) at the *sn*-1 and *sn*-3 positions and stearic acid (Item No. 10011298) at the *sn*-2 position. It has been found in soybean oil as well as transitional and mature human milk.^{1,2}

References

1. Gao, B., Luo, Y., Lu, W., *et al.* Triacylglycerol compositions of sunflower, corn and soybean oils examined with supercritical CO₂ ultra-performance convergence chromatography combined with quadrupole time-of-flight mass spectrometry. *Food Chem.* **218**, 569-574 (2017).
2. Tu, A., Ma, Q., Bai, H., *et al.* A comparative study of triacylglycerol composition in Chinese human milk within different lactation stages and imported infant formula by SFC coupled with Q-TOF-MS. *Food Chem.* **221**, 555-567 (2017).

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, 02/22/2024

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