

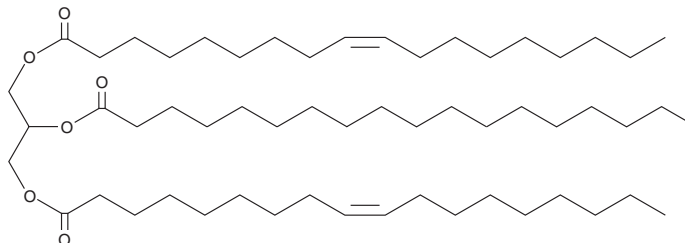
PRODUCT INFORMATION



1,3-Dioleoyl-2-Stearoyl Glycerol

Item No. 26889

CAS Registry No.: 2410-29-9
Formal Name: 9Z-octadecenoic acid, 1,1'-[2-[(1-oxooctadecyl)oxy]-1,3-propanediyl] ester
Synonyms: 1,3-Olein-2-Stearin, TG(18:1/18:0/18:1)
MF: C₅₇H₁₀₆O₆
FW: 887.5
Purity: ≥98%
Supplied as: A solid
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-Dioleoyl-2-stearoyl glycerol is supplied as a solid. A stock solution may be made by dissolving the 1,3-dioleoyl-2-stearoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Dioleoyl-2-stearoyl glycerol is slightly soluble in chloroform and methanol.

Description

1,3-Dioleoyl-2-stearoyl glycerol is a triacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 and *sn*-3 positions and stearic acid (Item No. 10011298) at the *sn*-2 position. It has been found in butterfat.¹

Reference

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).

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, 12/21/2022

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