

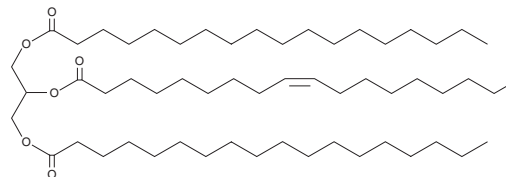
PRODUCT INFORMATION



1,3-Distearoyl-2-Oleoyl Glycerol

Item No. 27163

CAS Registry No.: 2846-04-0
Formal Name: (9Z)-9-Octadecenoic acid, 2-[(1-oxooctadecyl)oxy]-1-[[[(1-oxooctadecyl)oxy)methyl]ethyl ester
Synonyms: TG(18:0/18:1/18:0), 1,3-Distearin-2-Olein
MF: C₅₇H₁₀₈O₆
FW: 889.5
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-Distearoyl-2-oleoyl glycerol is supplied as a solid. A stock solution may be made by dissolving the 1,3-distearoyl-2-oleoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Distearoyl-2-oleoyl glycerol is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

1,3-Distearoyl-2-oleoyl glycerol is a triacylglycerol that contains stearic acid (Item No. 10011298) at the *sn*-1 and *sn*-3 positions and oleic acid (Item Nos. 90260 | 24659) at the *sn*-2 position. It has been found in Kpangnana, Chinese tallow, and cocoa butters.¹ 1,3-Distearoyl-2-oleoyl glycerol increases the phagocytic function of reticulo-endothelial cells in mice when administered at doses of 5 and 10 mg/animal.²

References

1. Ghazani, S.M. and Marangoni, A.G. The triclinic polymorphism of cocoa butter is dictated by its major molecular species, 1-palmitoyl, 2-oleoyl, 3-stearoyl glycerol (POS). *Cryst. Growth Des.* **19**, 90-97 (2019).
2. Stuart, A.E. and Cooper, G.N. Stimulation of the reticulo-endothelial phagocytic function by glyceryl tricaprinate and 2-oleodistearin. *Exp. Mol. Pathol.* **2(3)**, 215-218 (1963).

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, 12/16/2022

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