

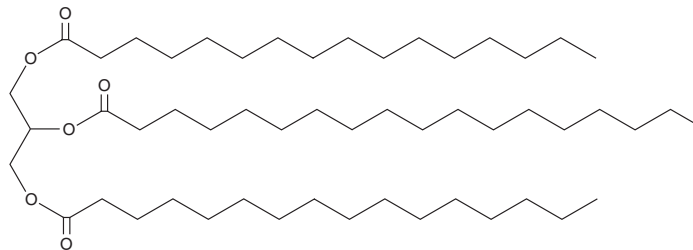
PRODUCT INFORMATION



1,3-Dipalmitoyl-2-Stearoyl Glycerol

Item No. 26989

CAS Registry No.: 2177-97-1
Formal Name: octadecanoic acid, 2-[[[1-oxohexadecyl)oxy]-1-[[[1-oxohexadecyl)oxy]methyl]ethyl ester
Synonyms: 1,3-Palmitin-2-Stearin, 16:0/18:0/16:0-TG, TG(16:0/18:0/16:0)
MF: C₅₃H₁₀₂O₆
FW: 835.4
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-Dipalmitoyl-2-stearoyl glycerol is supplied as a solid. A stock solution may be made by dissolving the 1,3-dipalmitoyl-2-stearoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Dipalmitoyl-2-stearoyl glycerol is soluble in the organic solvent dimethyl formamide at a concentration of approximately 10 mg/ml.

Description

1,3-Dipalmitoyl-2-stearoyl glycerol is a triacylglycerol that contains palmitic acid (Item No. 10006627) at the *sn*-1 and *sn*-3 positions and stearic acid (Item No. 10011298) at the *sn*-2 position. 1,3-Dipalmitoyl-2-stearoyl glycerol levels decrease in rats with reduced food intake-induced energy depletion compared to fed rats.¹

Reference

1. Chen, Z.-Y. and Cunnane, S.C. Short-term energy deficit causes net accumulation of linoleoyl-enriched triacylglycerols in rat liver. *FEBS Lett.* **280(2)**, 393-396 (1991).

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, 11/16/2022

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