

PRODUCT INFORMATION



1,2-Distearoyl-3-Palmitoyl-*rac*-glycerol

Item No. 28551

CAS Registry No.: 2177-99-3
Formal Name: octadecanoic acid,
1,1'-[1-[[[(1-oxohexadecyl)oxy]
methyl]-1,2-ethanediyl] ester

Synonyms: 1,2-Stearin-3-Palmitin,
TG(18:0/18:0/16:0)

MF: C₅₅H₁₀₆O₆

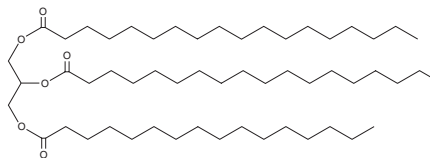
FW: 863.4

Purity: ≥95%

Supplied as: A crystalline 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

1,2-Distearoyl-3-palmitoyl-*rac*-glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-distearoyl-3-palmitoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1,2-Distearoyl-3-palmitoyl-*rac*-glycerol is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

1,2-Distearoyl-3-palmitoyl-*rac*-glycerol is a triacylglycerol that contains stearic acid (Item No. 10011298) at the *sn*-1 and *sn*-2 positions and palmitic acid (Item No. 10006627) at the *sn*-3 position. It has been found in cocoa butter, lard, and cod liver oil.^{1,2}

References

1. Buchgraber, M., Ulbeth, F., and Anklam, E. Comparison of HPLC and GLC techniques for the determination of the triglyceride profile of cocoa butter. *J. Agric. Food Chem.* **48(8)**, 3359-3363 (2000).
2. Rohman, A., Triyana, K., Sismindari, S., *et al.* Differentiation of lard and other animal fats based on triacylglycerols composition and principal component analysis. *Int. Food Res. J.* **19(2)**, 475-479 (2012).

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, 12/08/2022

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