

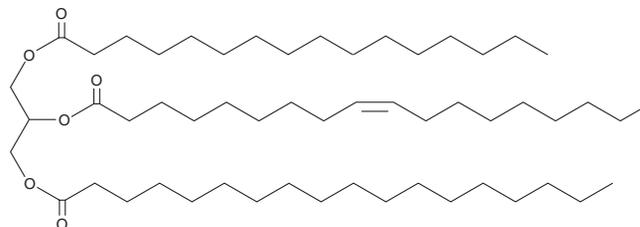
PRODUCT INFORMATION



1-Palmitoyl-2-Oleoyl-3-Stearoyl-*rac*-glycerol

Item No. 27067

CAS Registry No.: 2190-27-4
Formal Name: 9Z-octadecenoic acid, 1-[[[(1-oxohexadecyl)oxy]methyl]-2-[[[(1-oxooctadecyl)oxy]ethyl ester
Synonyms: 1-Palmitin-2-Olein-3-Stearin, TG(16:0/18:1/18:0)
MF: C₅₅H₁₀₄O₆
FW: 861.4
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-Palmitoyl-2-oleoyl-3-stearoyl-*rac*-glycerol is supplied as a solid. A stock solution may be made by dissolving the 1-palmitoyl-2-oleoyl-3-stearoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1-Palmitoyl-2-oleoyl-3-stearoyl-*rac*-glycerol is slightly soluble in chloroform and methanol.

Description

1-Palmitoyl-2-oleoyl-3-stearoyl-*rac*-glycerol is a triacylglycerol that contains palmitic acid (Item No. 10006627) at the *sn*-1 position, oleic acid (Item Nos. 90260 | 24659) at the *sn*-2 position, and stearic acid (Item No. 10011298) at the *sn*-3 position. It is one of the primary triacylglycerols found in cocoa butter.^{1,2}

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(1)**, 90-97 (2019).
2. Bootello, M.A., Hartel, R.W., Garcés, R., *et al.* Evaluation of high oleic-high stearic sunflower hard stearins for cocoa butter equivalent formulation. *Food Chem.* **134(3)**, 1409-1417 (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, 12/01/2022

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