

PRODUCT INFORMATION

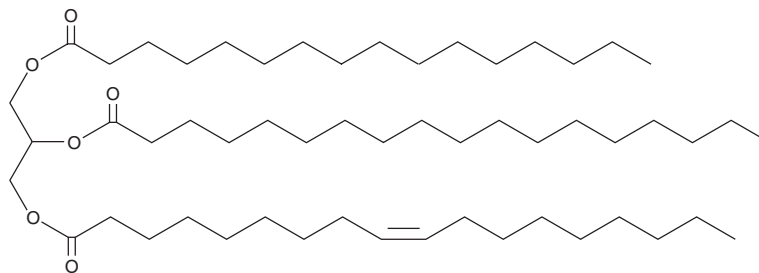


1-Palmitoyl-2-Stearoyl-3-Oleoyl-*rac*-glycerol

Item No. 27068

CAS Registry No.: 2190-28-5
Formal Name: 9Z-octadecenoic acid, 3-[(1-oxohexadecyl)oxy]-2-[(1-oxooctadecyl)oxy]propyl ester
Synonyms: 1-Palmitin-2-Stearin-3-Olein, TG(16:0/18:0/18:1)

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-stearoyl-3-oleoyl-*rac*-glycerol is supplied as a solid. A stock solution may be made by dissolving the 1-palmitoyl-2-stearoyl-3-oleoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1-Palmitoyl-2-stearoyl-3-oleoyl-*rac*-glycerol is slightly soluble in methanol and chloroform.

Description

1-Palmitoyl-2-stearoyl-3-oleoyl-*rac*-glycerol is a triacylglycerol that contains palmitic (Item No. 10006627), stearic (Item No. 10011298), and oleic acid (Item Nos. 90260 | 24659) at the *sn*-1, *sn*-2, and *sn*-3 positions, respectively. It has been detected in RAW 264.7 cells by neutral loss MS.¹ Increased serum levels of 1-palmitoyl-2-stearoyl-3-oleoyl-*rac*-glycerol are a potential biomarker for non-alcoholic fatty liver disease (NAFLD).²

References

1. Murphy, R.C., James, P.F., McAnoy, A.M., *et al.* Detection of the abundance of diacylglycerol and triacylglycerol molecular species in cells using neutral loss mass spectrometry. *Anal. Biochem.* **366**(1), 59-70 (2007).
2. Orešič, M., Hyötyläinen, T., Kotronen, A., *et al.* Prediction of non-alcoholic fatty-liver disease and liver fat content by serum molecular lipids. *Diabetologia* **56**(10), 2266-2274 (2013).

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, 12/20/2022

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