

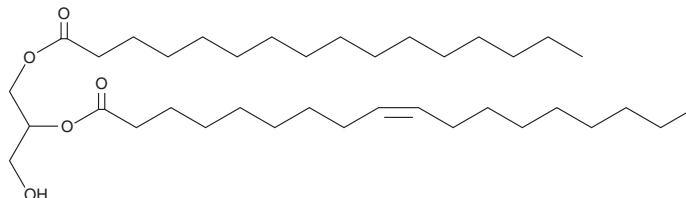
PRODUCT INFORMATION



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

Item No. 27164

CAS Registry No.: 3123-73-7
Formal Name: 9Z-octadecenoic acid, 1-(hydroxymethyl)-2-[(1-oxohexadecyl)oxy]ethyl ester
Synonyms: 16:0/18:1-DG, DG(16:0/18:1/0:0), 1-Palmitin-2-Olein
MF: C₃₇H₇₀O₅
FW: 595.0
Purity: ≥95%
Supplied as: An oil
Storage: -80°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-*rac*-glycerol is supplied as an oil. A stock solution may be made by dissolving the 1-palmitoyl-2-oleoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1-Palmitoyl-2-oleoyl-*rac*-glycerol is slightly soluble in chloroform and ethyl acetate.

Description

1-Palmitoyl-2-oleoyl-*rac*-glycerol is a diacylglycerol containing palmitic acid (Item No. 10006627) at the *sn*-1 position and oleic acid (Item Nos. 90260 | 24659) at the *sn*-2 position. It has been found in crude sunflower lecithin.¹ 1-Palmitoyl-2-oleoyl-*rac*-glycerol levels increase in RINm5F rat insulinoma cells treated with oleate, elaidate, or *trans*-vacenate.²

References

- Goñi, M.L., Pacheco, C., Constenla, D.T., *et al.* Solvent-free enzymatic hydrolysis of non-polar lipids in crude sunflower lecithin using phospholipase A₁ (Lecitase® Ultra). *Biocatal. Biotransfor.* **36(5)**, 341-351 (2017).
- Sarnyai, F., Donkó, M.B., Mátyási, J., *et al.* Cellular toxicity of dietary trans fatty acids and its correlation with ceramide and diglyceride accumulation. *Food Chem. Toxicol.* **124**, 324-335 (2019).

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, 01/21/2026

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