

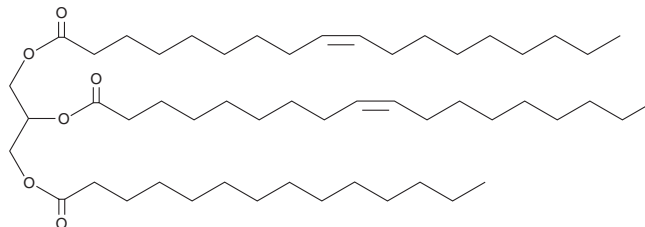
PRODUCT INFORMATION



1,2-Dioleoyl-3-Myristoyl-*rac*-glycerol

Item No. 27291

CAS Registry No.: 35804-98-9
Formal Name: 9Z-octadecenoic acid, 1,1'-[1-[[[1-oxotetradecyl]oxy]methyl]-1,2-ethanediyl] ester
Synonyms: 1,2-Olein-3-Myristin, TG(18:1/18:1/14:0)
MF: C₅₃H₉₈O₆
FW: 831.3
Purity: ≥98%
Supplied as: A 10 mg/ml solution in methyl acetate
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-Dioleoyl-3-myristoyl-*rac*-glycerol is supplied as a 10 mg/ml solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as chloroform and methanol purged with an inert gas can be used. 1,2-Dioleoyl-3-myristoyl-*rac*-glycerol is slightly soluble in these solvents.

Description

1,2-Dioleoyl-3-myristoyl-*rac*-glycerol is a triacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 and *sn*-2 positions and myristic acid (Item No. 13351) at the *sn*-3 position. It has been found in the fat body of male *B. lapidarius* bumblebees and in *R. erythropolis* bacteria utilizing phenol or succinic acid as a carbon source.^{1,2}

References

1. Cvačka, J., Hovorka, O., Jiroš, P., *et al.* Analysis of triacylglycerols in fat body of bumblebees by chromatographic methods. *J. Chromatogr. A*. **1101(1-2)**, 226-237 (2006).
2. Řezanka, T., Schreiberová, O., Krulíková, T., *et al.* RP-HPLC/MS-APCI analysis of odd-chain TAGs from *Rhodococcus erythropolis* including some regioisomers. *Chem. Phys. Lipids* **163(4-5)**, 373-380 (2010).

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, 12/03/2025

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