

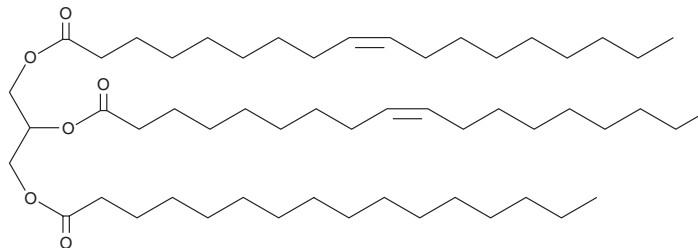
# PRODUCT INFORMATION



## 1,2-Dioleoyl-3-Palmitoyl-*rac*-glycerol

Item No. 27070

**CAS Registry No.:** 2190-30-9  
**Formal Name:** 9Z-octadecenoic acid  
1,1'-[1-[[[1-oxohexadecyl)oxy]  
methyl]-1,2-ethanediyl] ester  
**Synonyms:** 1,2-Olein-3-Palmitin, OOP,  
TG(18:1(9Z)/18:1(9Z)/16:0)  
**MF:** C<sub>55</sub>H<sub>102</sub>O<sub>6</sub>  
**FW:** 859.4  
**Purity:** ≥90%  
**Supplied as:** A solution in methyl acetate  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

1,2-Dioleoyl-3-palmitoyl-*rac*-glycerol is supplied as a 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. 1,2-Dioleoyl-3-palmitoyl-*rac*-glycerol is slightly soluble in chloroform and methanol.

### Description

1,2-Dioleoyl-3-palmitoyl-*rac*-glycerol is a triacylglycerol that contains oleic acid (Item No. 90260) at the *sn*-1 and *sn*-2 positions and palmitic acid (Item No. 10006627) at the *sn*-3 position. It has been found in *Leccino*, *Caroleo*, *Dritta*, and *Coratina* olive oils.<sup>1</sup>

### Reference

1. Ranalli, A., Lucera, L., Contento, S., *et al.* Bioactive constituents, flavors and aromas of virgin oils obtained by processing olives with a natural enzyme extract. *Eur. J. Lipid Sci. Technol.* **106(3)**, 187-197 (2004).

#### 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, 10/24/2022

#### 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