

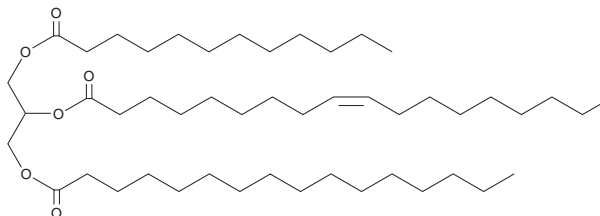
# PRODUCT INFORMATION



## 1-Lauroyl-2-Oleoyl-3-Palmitoyl-*rac*-glycerol

Item No. 28549

**CAS Registry No.:** 51604-51-4  
**Formal Name:** 9-octadecenoic acid (9Z)-1-[[[1-oxododecyl)oxy)methyl]-2-[[[1-oxohexadecyl)oxy]ethyl ester  
**Synonyms:** TG(12:0/18:1/16:0), 1-Laurin-2-Olein-3-Palmitin  
**MF:** C<sub>49</sub>H<sub>92</sub>O<sub>6</sub>  
**FW:** 777.3  
**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-Lauroyl-2-oleoyl-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-Lauroyl-2-oleoyl-3-palmitoyl-*rac*-glycerol is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

1-Lauroyl-2-oleoyl-3-palmitoyl-*rac*-glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 1-lauroyl-2-oleoyl-3-palmitoyl-*rac*-glycerol should be diluted with the aqueous buffer of choice. 1-Lauroyl-2-oleoyl-3-palmitoyl-*rac*-glycerol has a solubility of 0.25 mg/ml in a 1:3 solution of chloroform:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

1-Lauroyl-2-oleoyl-3-palmitoyl-*rac*-glycerol is a triacylglycerol that contains lauric acid (Item No. 10006626), oleic acid (Item Nos. 90260 | 24659), and palmitic acid (Item No. 10006627) at the *sn*-1, *sn*-2, and *sn*-3 positions, respectively. It has been found in date seed oil.<sup>1</sup>

### Reference

1. Holcapek, M., Lísá, M., Jandera, P., *et al.* Quantitation of triacylglycerols in plant oils using HPLC with APCI-MS, evaporative light-scattering, and UV detection. *J. Sep. Sci.* **28(12)**, 1315-1333 (2005).

#### 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, 09/09/2025

#### 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