

# 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:  $C_{49}H_{92}O_6$

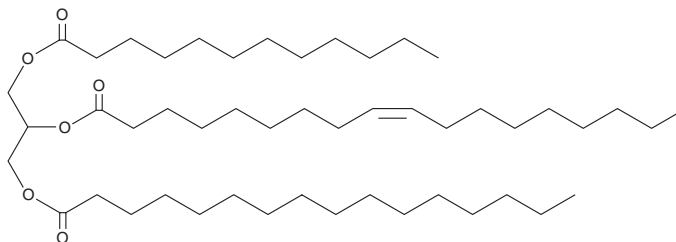
FW: 777.3

Purity:  $\geq 98\%$

Supplied as: A solid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  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 solid. A stock solution may be made by dissolving the 1-lauroyl-2-oleoyl-3-palmitoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 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, 1-lauroyl-2-oleoyl-3-palmitoyl-*rac*-glycerol should first be dissolved in chloroform and then diluted with the aqueous buffer of choice. 1-Lauroyl-2-oleoyl-3-palmitoyl-*rac*-glycerol has a solubility of approximately 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, 11/15/2022

#### 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