

# PRODUCT INFORMATION



## 1-Linoleoyl-3- $\alpha$ -Linolenoyl-*rac*-glycerol

Item No. 26829

**CAS Registry No.:** 126374-41-2  
**Formal Name:** 9Z,12Z,15Z-octadecatrienoic acid, 2-hydroxy-3-[[[(9Z,12Z)-1-oxo-9,12-octadecadien-1-yl]oxy]propyl ester  
**Synonyms:** DG(18:2/0:0/18:3), 1-Linolein-3- $\alpha$ -Linolenin, 1-Linolein-3-Linolenin

**MF:** C<sub>39</sub>H<sub>66</sub>O<sub>5</sub>

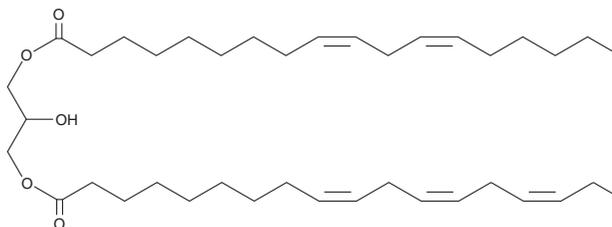
**FW:** 614.9

**Purity:**  $\geq$ 98%

**Supplied as:** A solution in ethanol

**Storage:** -20°C

**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

1-Linoleoyl-3- $\alpha$ -linolenoyl-*rac*-glycerol is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as methanol purged with an inert gas can be used. 1-Linoleoyl-3- $\alpha$ -linolenoyl-*rac*-glycerol is soluble in methanol.

### Description

1-Linoleoyl-3- $\alpha$ -linolenoyl-*rac*-glycerol is a diacylglycerol that contains linoleic acid (Item Nos. 90150 | 90150.1 | 21909) at the *sn*-1 position and  $\alpha$ -linolenic acid (Item Nos. 90210 | 21910) at the *sn*-3 position. It has been found in olive oil subjected to lipase-catalyzed glycerolysis with immobilized lipase B from *C. antarctica*.<sup>1</sup>

### Reference

1. Singh, A.K. and Mukhopadhyay, M. Olive oil glycerolysis with the immobilized lipase *Candida antarctica* in a solvent free system. *Grasas Y Aceites* **63(2)**, 202-208 (2012).

#### 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, 11/08/2024

#### 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