

PRODUCT INFORMATION

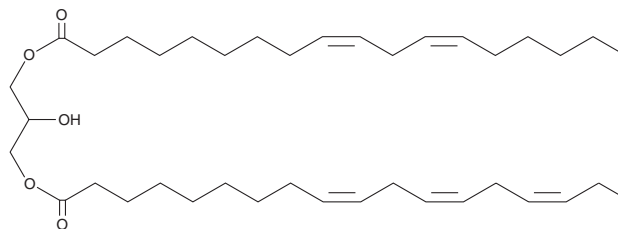


1-Linoleoyl-3- α -Linolenoyl-*rac*-glycerol

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- α -Linolenin, 1-Linolein-3-Linolenin

MF: C₃₉H₆₆O₅
FW: 614.9
Purity: \geq 98%
Supplied as: A neat oil
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- α -linolenoyl-*rac*-glycerol is supplied as a neat oil. A stock solution may be made by dissolving the 1-linoleoyl-3- α -linolenoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1-Linoleoyl-3- α -linolenoyl-*rac*-glycerol is soluble in methanol.

Description

1-Linoleoyl-3- α -linolenoyl-*rac*-glycerol is a diacylglycerol that contains linoleic acid (Item Nos. 90150 | 90150.1 | 21909) at the *sn*-1 position and α -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*.¹

References

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, 11/11/2022

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