

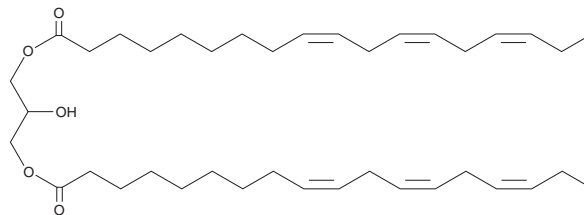
PRODUCT INFORMATION



1,3-Di- α -Linolenoyl Glycerol

Item No. 26940

CAS Registry No.: 98242-77-4
Formal Name: (9Z,9'Z,12Z,12'Z,15Z,15'Z)-9,12,15-octadecatrienoic acid, 2-hydroxy-1,3-propanediyl ester
Synonyms: DG(18:3/0:0/18:3), 1,3-Dilinolenin
MF: C₃₉H₆₄O₅
FW: 612.9
Purity: $\geq 95\%$
Supplied as: An oil
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

1,3-Di- α -linolenoyl glycerol is supplied as an oil. A stock solution may be made by dissolving the 1,3-di- α -linolenoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Di- α -linolenoyl glycerol is slightly soluble in chloroform and methanol.

Description

1,3-Di- α -linolenoyl glycerol is a diacylglycerol that contains α -linolenic acid (Item Nos. 90210 | 21910) at the *sn*-1 and *sn*-3 positions. It has been found in the roots of *A. sinensis* and the aerial part of *H. halodendron*.^{1,2}

References

1. Deng, S., Chen, S.-N., Yao, P., *et al.* Serotonergic activity-guided phytochemical investigation of the roots of *Angelica sinensis*. *J. Nat. Prod.* **69**(4), 536-541 (2006).
2. Kozhamkulova, Z.A., Radwan, M.M., Zhusupova, G.E., *et al.* Halimodendrin I, a new acylated triterpene glycoside from *Halimodendron halodendron* (Fabaceae). *Phytochem. Lett.* **4**(3), 323-327 (2011).

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, 01/26/2023

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