

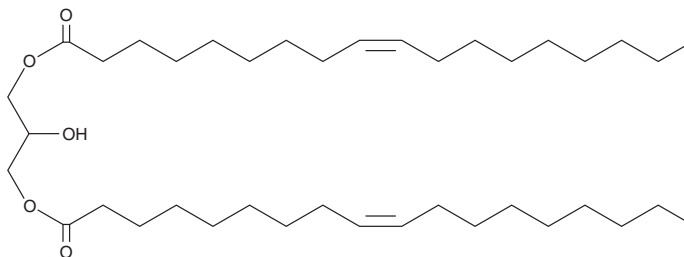
PRODUCT INFORMATION



1,3-Dioleoyl Glycerol

Item No. 26896

CAS Registry No.: 2465-32-9
Formal Name: 9Z-octadecenoic acid, 1,1'-(2-hydroxy-1,3-propanediyl) ester
Synonyms: DG(18:1/0:0/18:1), 1,3-Diolein
MF: C₃₉H₇₂O₅
FW: 621.0
Purity: ≥98%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

1,3-Dioleoyl glycerol is supplied as a liquid. A stock solution may be made by dissolving the 1,3-dioleoyl glycerol in the solvent of choice. 1,3-Dioleoyl glycerol is soluble in the organic solvent chloroform, which should be purged with an inert gas.

Description

1,3-Dioleoyl glycerol is a diacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 and *sn*-3 positions. It activates protein kinase C (PKC) by approximately 30% when used at a concentration of 50 μM.¹

Reference

1. Sánchez-Piñera, P., Micol, V., Corbalán-García, S., *et al.* A comparative study of the activation of protein kinase C α by different diacylglycerol isomers. *Biochem J.* **337**(3), 387-395 (1999).

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, 06/05/2020

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