

# PRODUCT INFORMATION



## 1,3-Dielaidoyl Glycerol

Item No. 26939

CAS Registry No.: 98168-52-6

Formal Name: (9E,9'E)-9-octadecenoic acid,  
1,1'-(2-hydroxy-1,3-propanediyl) ester

Synonyms: DG(18:1(9E)/0:0/18:1(9E)),  
DG(18:1/0:0/18:1), 1,3-Dielaidin

MF:  $C_{39}H_{72}O_5$

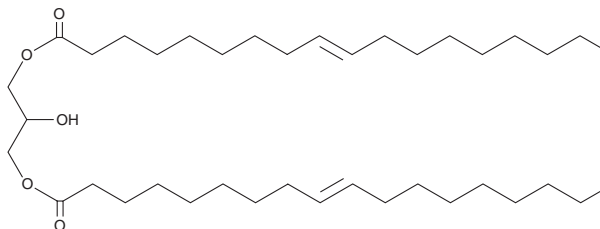
FW: 621.0

Purity:  $\geq 98\%$

Supplied as: An oil

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

1,3-Dielaidoyl glycerol is supplied as an oil. A stock solution may be made by dissolving the 1,3-dielaidoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Dielaidoyl glycerol is slightly soluble in organic solvents such as chloroform and ethyl acetate.

### Description

1,3-Dielaidoyl glycerol is a diacylglycerol that contains elaidic acid (Item No. 90250) at the *sn*-1 and *sn*-3 positions.

#### 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, 11/10/2022

#### 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