

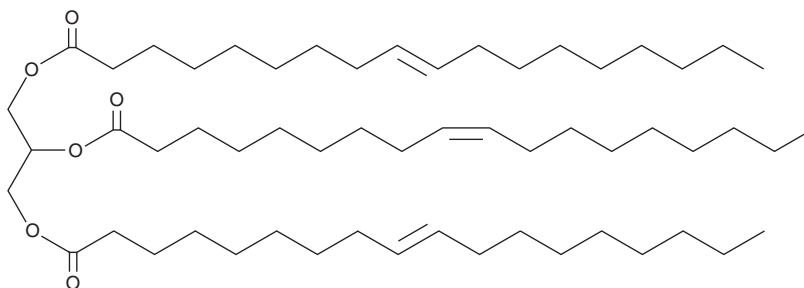
PRODUCT INFORMATION



1,3-Dielaidoyl-2-Oleoyl Glycerol

Item No. 29236

CAS Registry No.: 946577-28-2
Formal Name: (9E,9'E)-9-octadecenoic acid, 2-[[[(9Z)-1-oxo-9-octadecen-1-yl]oxy]-1,3-propanediyl ester
Synonyms: 1,3-Elaidin-2-Olein, TG(18:1/18:1/18:1)
MF: C₅₇H₁₀₄O₆
FW: 885.4
Purity: ≥95%
Supplied as: A neat 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-Dielaidoyl-2-oleoyl glycerol is supplied as a neat oil. A stock solution may be made by dissolving the 1,3-dielaidoyl-2-oleoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Dielaidoyl-2-oleoyl glycerol is miscible in organic solvents such as DMSO and dimethyl formamide.

Description

1,3-Dielaidoyl-2-oleoyl glycerol is a triacylglycerol that contains elaidic acid (Item No. 90250) at the *sn*-1 and *sn*-3 positions and oleic acid (Item Nos. 90260 | 24659) at the *sn*-2 position.

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, 03/01/2023

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