

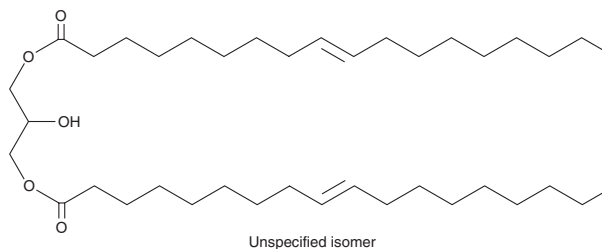
PRODUCT INFORMATION



Dielaidin

Item No. 26979

CAS Registry No.: 372516-91-1
Formal Name: (9E,9'E)-9-octadecenoic acid, diester with 1,2,3-propanetriol
MF: C₃₉H₇₂O₅
FW: 621.0
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Dielaidin is supplied as a solid. A stock solution may be made by dissolving the dielaidin in the solvent of choice, which should be purged with an inert gas. Dielaidin is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of dielaidin in these solvents is approximately 10 mg/ml.

Dielaidin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, dielaidin should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Dielaidin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Dielaidin is a diacylglycerol that contains the monounsaturated 18-carbon fatty acid elaidic acid (Item No. 90250), the *trans* isomer of oleic acid (Item Nos. 90260 | 24659), at two 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, 11/30/2022

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