

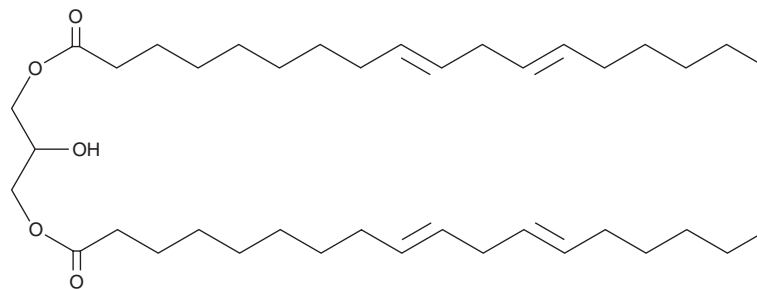
PRODUCT INFORMATION



Dilinoelaidin

Item No. 26980

CAS Registry No.: 372516-93-3
Formal Name: (9E,9'E,12E,12'E)-9,12-octadecadienoic acid, diester with 1,2,3-propanetriol
MF: C₃₉H₆₈O₅
FW: 617.0
Purity: ≥98%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



Unspecified isomer

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

Laboratory Procedures

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

Dilinoelaidin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, dilinoelaidin should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Dilinoelaidin 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

Dilinoelaidin is a diacylglycerol that contains the polyunsaturated 18-carbon fatty acid linoelaidic acid (Item No. 90160) 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, 08/08/2019

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