

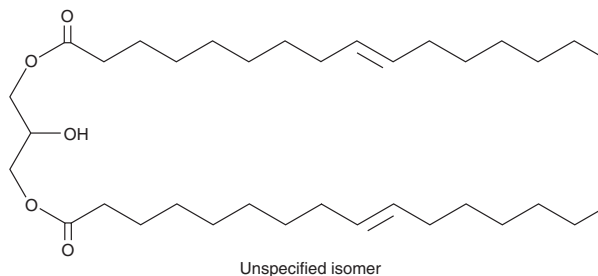
PRODUCT INFORMATION



Dipalmitelaidin

Item No. 26978

CAS Registry No.: 372516-90-0
Formal Name: (9E,9'E)-9-hexadecenoic acid, diester with 1,2,3-propanetriol
MF: C₃₅H₆₄O₅
FW: 564.9
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

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

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

Dipalmitelaidin is a diacylglycerol that contains the monounsaturated 16-carbon fatty acid palmitelaidic acid (Item No. 9001798) 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, 12/21/2022

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