

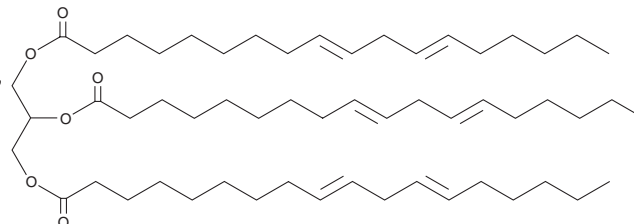
PRODUCT INFORMATION



1,2,3-Trilinoelaidoyl Glycerol

Item No. 26948

CAS Registry No.: 5188-25-0
Formal Name: (9E,9'E,9''E,12E,12'E,12''E)-9,12-octadecadienoic acid, 1,1',1''-(1,2,3-propanetriyl) ester
Synonyms: TG(18:2(9E,12E)/18:2(9E,12E)/18:2(9E,12E)), Trilinoelaidin
MF: C₅₇H₉₈O₆
FW: 879.4
Purity: ≥98%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

1,2,3-Trilinoelaidoyl glycerol is supplied as a liquid. A stock solution may be made by dissolving the 1,2,3-trilinoelaidoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Trilinoelaidoyl glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 1,2,3-trilinoelaidoyl glycerol in these solvents is approximately 10 mg/ml.

1,2,3-Trilinoelaidoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,2,3-trilinoelaidoyl glycerol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 1,2,3-Trilinoelaidoyl glycerol 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

1,2,3-Trilinoelaidoyl glycerol is a triacylglycerol that contains linoelaidic acid (Item No. 90160) at the *sn*-1, *sn*-2, and *sn*-3 positions. Dietary administration of 1,2,3-trilinoelaidoyl glycerol (0.6-6.3% w/w) lowers serum levels of thromboxane B₂ (TXB₂; Item No. 19030), prostaglandin F₂ (PGF₂), and PGE in rats.¹

Reference

1. Bruckner, G., Goswami, S., and Kinsella, J.E. Dietary trilinoelaidate: Effects on organ fatty acid composition, prostanoid biosynthesis and platelet function in rats. *J. Nutr.* **114**(1), 58-67 (1984).

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, 06/05/2020

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