

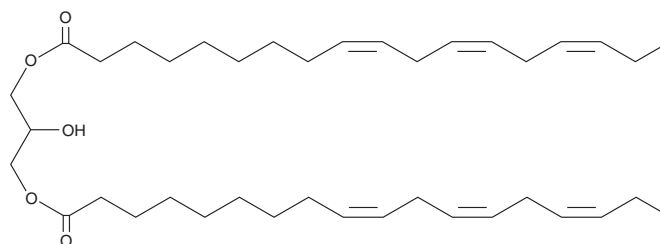
PRODUCT INFORMATION



Di- α -linolenin

Item No. 27013

CAS Registry No.: 195711-48-9
Formal Name: (9Z,9'Z,12Z,12'Z,15Z,15'Z)-9,12,15-octadecatrienoic acid, diester with 1,2,3-propanetriol
Synonym: Dilinolenin
MF: C₃₉H₆₄O₅
FW: 612.9
Purity: \geq 95%
Supplied as: A liquid
Storage: -20°C
Stability: \geq 2 years



Unspecified isomer

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

Laboratory Procedures

Di- α -linolenin is supplied as a liquid. A stock solution may be made by dissolving the di- α -linolenin in the solvent of choice, which should be purged with an inert gas. Di- α -linolenin is soluble in organic solvents such as chloroform, ethanol, DMSO, dimethyl formamide (DMF). The solubility of di- α -linolenin in chloroform is approximately 10 mg/ml and approximately 30 mg/ml in ethanol, DMSO, and DMF.

Description

Di- α -linolenin is a diacylglycerol that contains the polyunsaturated 18-carbon fatty acid α -linolenic acid (Item Nos. 90210 | 21910) 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, 01/16/2020

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