

PRODUCT INFORMATION



1,2,3-Tri- γ -Linolenoyl Glycerol

Item No. 10009825

CAS Registry No.: 60756-74-3
Formal Name: (6Z,6'Z,6''Z,9Z,9'Z,9''Z,12Z,12'Z,12''Z)-6,9,12-octadecatrienoic acid, 1,1,1''-(1,2,3-propanetriyl) ester
Synonyms: γ -Linolenic Acid Triglyceride, γ -Linolenin, TG(18:3/18:3/18:3), Tri- γ -linolenoate

MF: C₅₇H₉₂O₆

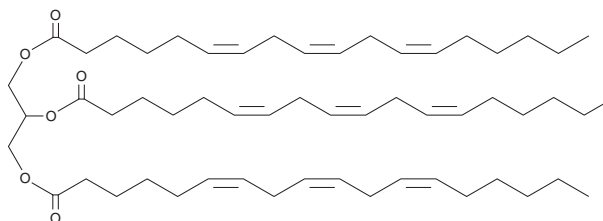
FW: 873.4

Purity: $\geq 98\%$

Supplied as: A solution in ethanol

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-Tri- γ -linolenoyl glycerol is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. 1,2,3-Tri- γ -linolenoyl glycerol is soluble in the organic solvent dimethyl formamide, which should be purged with an inert gas, at a concentration of approximately 10 mg/ml.

Description

1,2,3-Tri- γ -linolenoyl glycerol is a polyunsaturated triacylglycerol with γ -linolenoyl (Item No. 90220) side chains attached at the C-1, C-2, and C-3 positions. It has been found in various plant species, including blackcurrant, redcurrant, and borage.¹

Reference

1. Lísá, M. and Holcapek, M. Triacylglycerols profiling in plant oils important in food industry, dietetics and cosmetics using high-performance liquid chromatography-atmospheric pressure chemical ionization mass spectrometry. *J. Chromatogr. A* **1198-1199**, 115-130 (2008).

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, 06/04/2020

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