

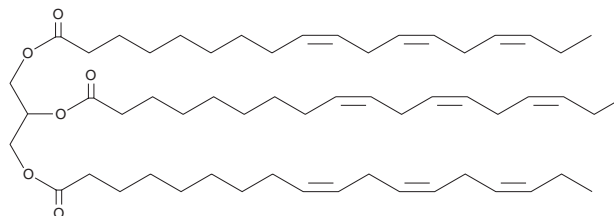
# PRODUCT INFORMATION



## 1,2,3-Tri- $\alpha$ -Linolenoyl Glycerol

Item No. 10009758

**CAS Registry No.:** 14465-68-0  
**Formal Name:** (9Z,9'Z,9''Z,12Z,12'Z,12''Z,15Z,15'Z,15''Z)-9,12,15-octadecatrienoic acid, 1,1',1''-(1,2,3-propanetriyl) ester  
**Synonyms:** Glycerol Trilinolenate, Glyceryl Trilinolenate, TG(18:3/18:3/18:3), Trilinolenin, Tri- $\alpha$ -linolenate  
**MF:** C<sub>57</sub>H<sub>92</sub>O<sub>6</sub>  
**FW:** 873.3  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:**  $\geq$ 2 years



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

### Laboratory Procedures

1,2,3-Tri- $\alpha$ -linolenoyl glycerol is supplied as a solution in ethanol. To change the solvent, simply evaporate the 1,2,3-tri- $\alpha$ -linolenoyl glycerol under a gentle stream of nitrogen and immediately add the solvent of choice. The solvent dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of 1,2,3-tri- $\alpha$ -linolenoyl glycerol in DMF is approximately 10 mg/ml.

1,2,3-Tri- $\alpha$ -linolenoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 1,2,3-tri- $\alpha$ -linolenoyl glycerol should be diluted with the aqueous buffer of choice. 1,2,3-Tri- $\alpha$ -linolenoyl glycerol has a solubility of approximately 500  $\mu$ g/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-Tri- $\alpha$ -linolenoyl glycerol is a polyunsaturated triacylglycerol with  $\alpha$ -linolenoyl (18:3) side-chains attached at the C-1, C-2, and C-3 positions that is useful in lipid research.

#### 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, 11/08/2022

#### 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](http://WWW.CAYMANCHEM.COM)