

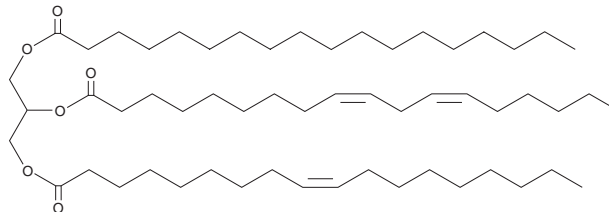
# PRODUCT INFORMATION



## 1-Stearoyl-2-Linoleoyl-3-Oleoyl-*rac*-glycerol

Item No. 28562

**CAS Registry No.:** 7162-26-7  
**Formal Name:** 9Z,12Z-octadecadienoic acid, 1-[[[(9Z)-1-oxo-9-octadecen-1-yl]oxy]methyl]-2-[(1-oxooctadecyl)oxy]ethyl ester  
**Synonyms:** 1-Stearin-2-Linolein-3-Olein, TG(18:0/18:2/18:1)  
**MF:** C<sub>57</sub>H<sub>104</sub>O<sub>6</sub>  
**FW:** 885.4  
**Purity:** ≥98%  
**Supplied as:** A solution in methyl acetate  
**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-Stearoyl-2-linoleoyl-3-oleoyl-*rac*-glycerol is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol and dimethyl formamide purged with an inert gas can be used. The solubility of 1-stearoyl-2-linoleoyl-3-oleoyl-*rac*-glycerol in these solvents is approximately 10 mg/ml.

For maximum solubility in aqueous buffers, evaporate the methyl acetate and dissolve in ethanol. The ethanolic solution of 1-stearoyl-2-linoleoyl-3-oleoyl-*rac*-glycerol should be diluted with the aqueous buffer of choice. 1-Stearoyl-2-linoleoyl-3-oleoyl-*rac*-glycerol has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method.

### Description

1-Stearoyl-2-linoleoyl-3-oleoyl-*rac*-glycerol is a triacylglycerol that contains stearic acid (Item No. 10011298), linoleic acid (Item Nos. 90150 | 90150.1 | 21909), and oleic acid (Item Nos. 90260 | 24659) at the *sn*-1, *sn*-2, and *sn*-3 positions, respectively. It has been found in poppy seed, maize, and evening primrose oils.<sup>1</sup>

### Reference

1. Mottram, H.R., Woodbury, S.E., and Evershed, R.P. Identification of triacylglycerol positional isomers present in vegetable oils by high performance liquid chromatography/atmospheric pressure chemical ionization mass spectrometry. *Rapid Commun. Mass Spectrom.* **11(12)**, 1240-1252 (1997).

#### 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, 12/02/2020

#### 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