

# PRODUCT INFORMATION



## 1,2-Di-13(Z)-Docosenoyl-3-Oleoyl-*rac*-glycerol

Item No. 27162

CAS Registry No.: 60003-02-3

Formal Name: (Z,Z,Z)-13-docosenoic acid, 1-[[[(1-oxo-9-octadecenyl)oxy]methyl]-1,2-ethanediyl ester

Synonyms: 1,2-13(Z)-Docosenoyl-3-Oleoyl-*rac*-glycerol, 1,2-Erucin(13Z)-3-Olein, TG(22:1(13Z)/22:1(13Z)/18:1(9Z))

MF:  $C_{65}H_{120}O_6$

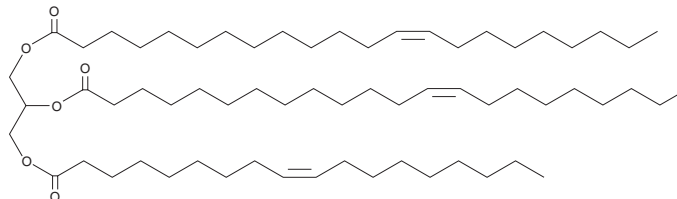
FW: 997.6

Purity:  $\geq 98\%$

Supplied as: A liquid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 2$  years



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

### Laboratory Procedures

1,2-Di-13(Z)-docosenoyl-3-oleoyl-*rac*-glycerol is supplied as a liquid. A stock solution may be made by dissolving the 1,2-di-13(Z)-docosenoyl-3-oleoyl-*rac*-glycerol in the solvent of choice. 1,2-Di-13(Z)-docosenoyl-3-oleoyl-*rac*-glycerol is soluble in organic solvents such as ethanol and dimethyl formamide, which should be purged with an inert gas. The solubility of 1,2-di-13(Z)-docosenoyl-3-oleoyl-*rac*-glycerol in these solvents is approximately 10 mg/ml.

1,2-Di-13(Z)-docosenoyl-3-oleoyl-*rac*-glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,2-di-13(Z)-docosenoyl-3-oleoyl-*rac*-glycerol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 1,2-Di-13(Z)-docosenoyl-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. We do not recommend storing the aqueous solution for more than one day.

### Description

1,2-Di-13(Z)-docosenoyl-3-oleoyl-*rac*-glycerol is a triacylglycerol that contains 13(Z)-docosenoic acid (Item No. 90175) at the *sn*-1 and *sn*-2 positions and oleic acid (Item No. 90260) at the *sn*-3 position.

#### 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, 02/19/2019

#### 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