

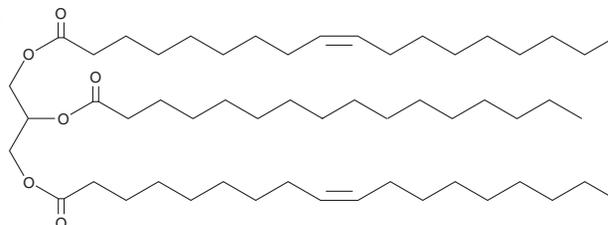
# PRODUCT INFORMATION



## 1,3-Dioleoyl-2-Palmitoyl Glycerol

Item No. 26843

**CAS Registry No.:** 1716-07-0  
**Formal Name:** 9(Z)-octadecenoic acid, 1,1'-[2-[(1-oxohexadecyl)oxy]-1,3-propanediyl] ester  
**Synonyms:** 1,3-Olein-2-Palmitin, TG(18:1/16:0/18:1)  
**MF:** C<sub>55</sub>H<sub>102</sub>O<sub>6</sub>  
**FW:** 859.4  
**Purity:** ≥98%  
**Supplied as:** A solution in hexane  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

1,3-Dioleoyl-2-palmitoyl glycerol is supplied as a solution in hexane. A stock solution may be made by dissolving the 1,3-dioleoyl-2-palmitoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Dioleoyl-2-palmitoyl glycerol is soluble in the organic solvent chloroform.

### Description

1,3-Dioleoyl-2-palmitoyl glycerol is a triacylglycerol with oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 and *sn*-3 positions and palmitic acid (Item No. 10006627) at the *sn*-2 position. It is one of the most common triacylglycerols found in human breast milk and levels are higher in colostrum than transitional and mature milk.<sup>1</sup> Human breast milk levels of 1,3-dioleoyl-2-palmitoyl glycerol are higher in Western women compared with Chinese women.

### Reference

1. Yuan, T.L., Qi, C., Dai, X., *et al.* Triacylglycerol composition of breast milk during different lactation stages. *J. Agric. Food Chem.* **67**(8), 2272-2278 (2019).

#### 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, 10/19/2022

#### 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