

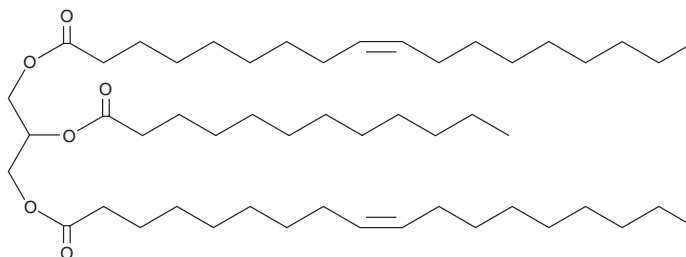
# PRODUCT INFORMATION



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

Item No. 27287

**CAS Registry No.:** 383189-71-7  
**Formal Name:** (9Z)-9-octadecenoic acid, 1,1'-[2-[(1-oxododecyl)oxy]-1,3-propanediyl] ester  
**Synonyms:** 1,3-Olein-2-Laurin, TG(18:1/12:0/18:1)  
**MF:** C<sub>51</sub>H<sub>94</sub>O<sub>6</sub>  
**FW:** 803.3  
**Purity:** ≥98%  
**Supplied as:** An oil  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

1,3-Dioleoyl-2-lauroyl glycerol is supplied as an oil. A stock solution may be made by dissolving the 1,3-dioleoyl-2-lauroyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Dioleoyl-2-lauroyl glycerol is slightly soluble in chloroform and methanol.

### Description

1,3-Dioleoyl-2-lauroyl glycerol is a triacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 and *sn*-3 positions and lauric acid (Item No. 10006626) at the *sn*-2 position. It has been found in human milk, butterfat, and palm kernel oil.<sup>1-3</sup>

### References

1. Giuffrida, F., Marmet, C., Tavazzi, I., et al. Quantification of 1,3-olein-2-palmitin (OPO) and palmitic acid in *sn*-2 position of triacylglycerols in human milk by liquid chromatography coupled with mass spectrometry. *Molecules* **24**(1), E22 (2018).
2. Small, D.M. The effects of glyceride structure on absorption and metabolism. *Annu. Rev. Nutr.* **11**, 413-434 (1991).
3. Kalo, P., Kemppinen, A., and Ollilainen, V. Determination of triacylglycerols in butterfat by normal-phase HPLC and electrospray-tandem mass spectrometry. *Lipids* **44**(2), 169-195 (2009).

#### 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, 12/22/2022

#### 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)