

PRODUCT INFORMATION



1,3-Dioleoyl-2-Decanoyl Glycerol

Item No. 27276

CAS Registry No.: 81246-69-7

Formal Name: 9Z-octadecenoic acid,
1,1'-[2-[(1-oxodecyl)oxy]-1,3-
propanediyl] ester

Synonyms: 1,3-Dioleoyl-2-Caproyl Glycerol,
1,3-Olein-2-Caprin, 1,3-Olein-2-
Decanoin, TG(18:1/10:0/18:1)

MF: $C_{49}H_{90}O_6$

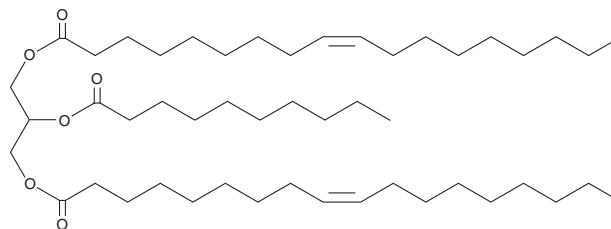
FW: 775.2

Purity: $\geq 98\%$

Supplied as: A liquid

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

Description

1,3-Dioleoyl-2-decanoyl glycerol is a triacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 and *sn*-3 positions and decanoic acid (Item No. 20838) at the *sn*-2 position. It has been found in transitional human milk.¹

Reference

1. Tu, A., Ma, Q., Bai, H., *et al.* A comparative study of triacylglycerol composition in Chinese human milk within different lactation stages and imported infant formula by SFC coupled with Q-TOF-MS. *Food Chem.* **221**, 555-567 (2017).

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, 06/08/2020

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