

PRODUCT INFORMATION



1,3-Dioleoyl-2-Octanoyl Glycerol

Item No. 26836

CAS Registry No.: 139979-77-4
Formal Name: 9Z-octadecenoic acid, 2-[(1-oxooctyl)oxy]-1,3-propanediyl ester

Synonyms: 1,3-Dioleoyl-2-Capryloyl Glycerol, OCO, 1,3-Olein-2-Caprylin, 1,3-Oleoyl-2-Capryloyl Glycerol, 1,3-Oleoyl-2-Octanoyl Glycerol, TG(18:1/8:0/18:1)

MF: $C_{47}H_{86}O_6$

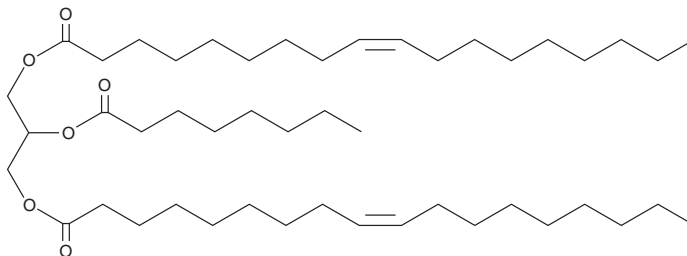
FW: 747.2

Purity: $\geq 98\%$

Supplied as: A liquid

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

Description

1,3-Dioleoyl-2-octanoyl glycerol is a triacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 and *sn*-3 positions and octanoic acid at the *sn*-2 position. It has been used to characterize the functionality of *R. oryzae* lipase for regiodistribution analysis of fats containing medium- and short-chain fatty acids.¹

Reference

1. Pérignon, M., Lecomte, J., Pina, M., et al. Evaluation of *Rhizopus oryzae* lipase for the determination of regiodistribution in triacylglycerols with medium chain fatty acids. *J. Am. Oil Chem. Soc.* **89**(1), 89-96 (2012).

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, 01/16/2026

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