

# PRODUCT INFORMATION



## 1,2-Dioctanoyl-3-Stearoyl-*rac*-glycerol

Item No. 27285

CAS Registry No.: 56149-05-4

Formal Name: octadecanoic acid, 2,3-bis[(1-oxooctyl)oxy]propyl ester

Synonyms: 1,2-Caprylin-3-Stearin,  
1,2-Octanoin-3-Stearin,  
TG(8:0/8:0/18:0)

MF:  $C_{37}H_{70}O_6$

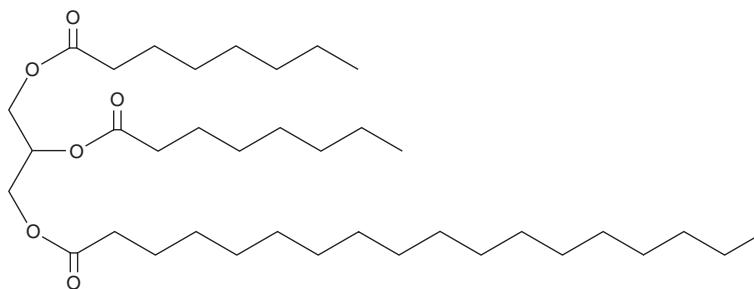
FW: 611.0

Purity:  $\geq 98\%$

Supplied as: An oil

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

Stability:  $\geq 4$  years



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

### Laboratory Procedures

1,2-Dioctanoyl-3-stearoyl-*rac*-glycerol is supplied as an oil. A stock solution may be made by dissolving the 1,2-dioctanoyl-3-stearoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1,2-Dioctanoyl-3-stearoyl-*rac*-glycerol is soluble in the organic solvent chloroform at a concentration of approximately 1 mg/ml.

### Description

1,2-Dioctanoyl-3-stearoyl-*rac*-glycerol is a triacylglycerol that contains octanoic acid at the *sn*-1 and *sn*-2 positions and stearic acid (Item No. 10011298) at the *sn*-3 position. It has been found in French butterfat.<sup>1</sup>

### Reference

1. Bornaz, S., Novak, G., and Parmentier, M. Seasonal and regional variation in triglyceride composition of French butterfat. *J. Am. Oil Chem. Soc.* **69**(11), 1131-1135 (1992).

#### 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