

PRODUCT INFORMATION



1,2-Dioleoyl-3-Docosanoyl-*rac*-glycerol

Item No. 28863

CAS Registry No.: 77145-66-5

Formal Name: docosanoic acid, 2,3-bis[[(9Z)-1-oxo-9-octadecen-1-yl]oxy]propyl ester

Synonyms: 1,2-Olein-3-Behenin,
TG(18:1/18:1/22:0)

MF: $C_{61}H_{114}O_6$

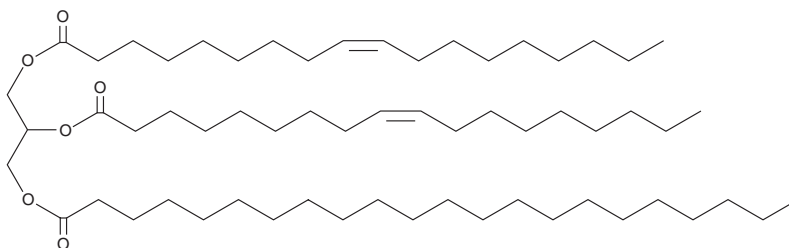
FW: 943.6

Purity: $\geq 95\%$

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥ 4 years



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

Laboratory Procedures

1,2-Dioleoyl-3-docosanoyl-*rac*-glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-dioleoyl-3-docosanoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1,2-Dioleoyl-3-docosanoyl-*rac*-glycerol is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml.

Description

1,2-Dioleoyl-3-docosanoyl-*rac*-glycerol is a triacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 and *sn*-2 positions and docosanoic acid (Item No. 9000338) at the *sn*-3 position. It has been found in cocoa butter and in cocoa butter equivalents (CBEs) containing shea stearin and palm mid fraction.¹

Reference

1. Bootello, M.A., Hartel, R.W., Garcés, R., *et al.* Evaluation of high oleic-high stearic sunflower hard stearins for cocoa butter equivalent formulation. *Food Chem.* **134**(3), 1409-1417 (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, 10/03/2022

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