

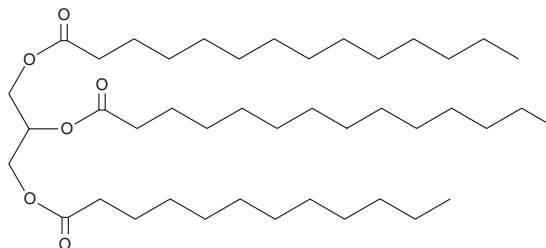
PRODUCT INFORMATION



1,2-Dimyristoyl-3-Lauroyl-*rac*-glycerol

Item No. 28861

CAS Registry No.: 60138-12-7
Formal Name: tetradecanoic acid, 1,1'-[1-[[[1-oxododecyl]oxy]methyl]-1,2-ethanediyl] ester
Synonyms: 1,2-Myristin-3-Laurin, TG(14:0/14:0/12:0)
MF: C₄₃H₈₂O₆
FW: 695.1
Purity: ≥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-Dimyristoyl-3-lauroyl-*rac*-glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-dimyristoyl-3-lauroyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1,2-Dimyristoyl-3-lauroyl-*rac*-glycerol is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml.

Description

1,2-Dimyristoyl-3-lauroyl-*rac*-glycerol is a triacylglycerol that contains myristic acid (Item No. 13351) at the *sn*-1 and *sn*-2 positions and lauric acid (Item No. 10006626) at the *sn*-3 position. It has been found in butterfat, aged cheese, and date seed oil.¹⁻³

References

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).
2. Nájera, A.I., Perea, S., Barcina, Y., *et al.* A reverse-phase high-performance liquid chromatographic method for analyzing complex mixtures of triglycerides. Application to the fat fraction of an aged cheese. *J. Am. Oil Chem. Soc.* **76(3)**, 399-407 (1999).
3. Holcapek, M., Lísá, M., Jandera, P., *et al.* Quantitation of triacylglycerols in plant oils using HPLC with APCI-MS, evaporative light-scattering, and UV detection. *J. Sep. Sci.* **28(12)**, 1315-1333 (2005).

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, 11/09/2022

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