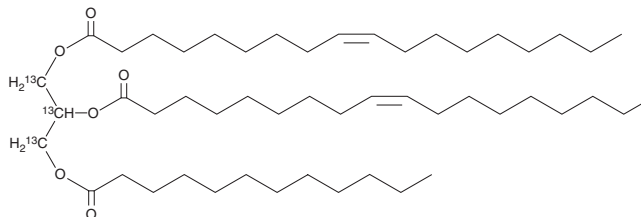


PRODUCT INFORMATION



1,2-Dioleoyl-3-Lauroyl-*rac*-glycerol-¹³C₃ Item No. 28755

CAS Registry No.: 2692624-36-3
Formal Name: 9Z-octadecenoic acid, 1,1'-[1-[[[1-oxododecyl]oxy]methyl-¹³C]-1,2-ethanediyl-1,2-¹³C₂] ester
Synonyms: 1,2-Olein-3-Laurin-¹³C₃, TG(18:1/18:1/12:0)-¹³C₃
MF: C₄₈[¹³C₃]H₉₄O₆
FW: 806.3
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

1,2-Dioleoyl-3-lauroyl-*rac*-glycerol-¹³C₃ is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as chloroform and methanol purged with an inert gas can be used. 1,2-Dioleoyl-3-lauroyl-*rac*-glycerol-¹³C₃ is slightly soluble in chloroform and methanol.

Description

1,2-Dioleoyl-3-lauroyl-*rac*-glycerol-¹³C₃ is intended for use as an internal standard for the quantification of 1,2-dioleoyl-3-lauroyl-*rac*-glycerol (Item No. 28548) by GC- or LC-MS. 1,2-Dioleoyl-3-lauroyl-*rac*-glycerol is a triacylglycerol that contains oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 and *sn*-2 positions and lauric acid (Item No. 10006626) at the *sn*-3 position. It has been found in date seed oil and the fat body of male *B. lapidarius* bumblebees.^{1,2}

References

- 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).
- Cvačka, J., Hovorka, O., Jiroš, P., *et al.* Analysis of triacylglycerols in fat body of bumblebees by chromatographic methods. *J. Chromatogr. A.* **1101(1-2)**, 226-237 (2006).

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, 12/02/2024

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