

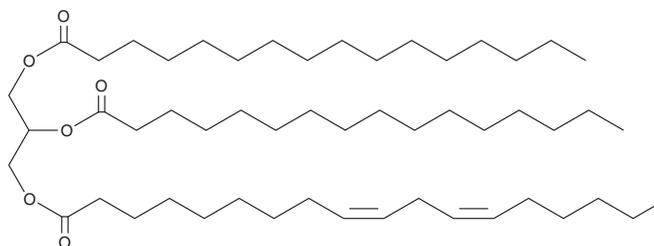
PRODUCT INFORMATION



1,2-Dipalmitoyl-3-Linoleoyl-*rac*-glycerol

Item No. 27160

CAS Registry No.: 2535-35-5
Formal Name: 9Z,12Z-octadecadienoic acid, 2,3-bis[(1-oxohexadecyl)oxy]propyl ester
Synonyms: 1,2-Palmitin-3-Linolein, TG(16:0/16:0/18:2)
MF: C₅₃H₉₈O₆
FW: 831.3
Purity: ≥98%
Supplied as: A 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-Dipalmitoyl-3-linoleoyl-*rac*-glycerol is supplied as a solid. A stock solution may be made by dissolving the 1,2-dipalmitoyl-3-linoleoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1,2-Dipalmitoyl-3-linoleoyl-*rac*-glycerol is slightly soluble in chloroform and methanol.

Description

1,2-Dipalmitoyl-3-linoleoyl-*rac*-glycerol is a triacylglycerol that contains palmitic acid (Item No. 10006627) at the *sn*-1 and *sn*-2 positions and linoleic acid (Item Nos. 90150 | 90150.1 | 21909) at the *sn*-3 position. It has been found in ostrich and emu oils and bovine milk fat.^{1,2}

References

1. Zhou, Y., Xue, Y., Chen, G.C., *et al.* Rapid separation and characterisation of triacylglycerols in ostrich oil by ultra performance liquid chromatography coupled with quadrupole time-of-flight mass spectrometry. *Food Chem.* **141**(3), 2098-2102 (2013).
2. Zhou, Q., Gao, B., Zhang, X., *et al.* Chemical profiling of triacylglycerols and diacylglycerols in cow milk fat by ultra-performance convergence chromatography combined with a quadrupole time-of-flight mass spectrometry. *Food Chem.* **143**, 199-204 (2014).

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, 12/15/2022

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