

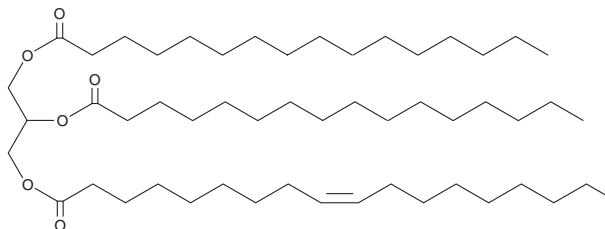
PRODUCT INFORMATION



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

Item No. 26878

CAS Registry No.: 1867-91-0
Formal Name: 9Z-octadecenoic acid, 2,3-bis[(1-oxohexadecyl)oxy]propyl ester
Synonyms: 1,2-Palmitin-3-Olein, TG(16:0/16:0/18:1)
MF: C₅₃H₁₀₀O₆
FW: 833.4
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-Dipalmitoyl-3-oleoyl-*rac*-glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-dipalmitoyl-3-oleoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1,2-Dipalmitoyl-3-oleoyl-*rac*-glycerol is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

1,2-Dipalmitoyl-3-oleoyl-*rac*-glycerol is a triacylglycerol that contains palmitic acid (Item No. 10006627) at the *sn*-1 and *sn*-2 positions and oleic acid (Item Nos. 90260 | 24659) at the *sn*-3 position. It has been found in *Apiaceae* seed oils, lard, and tallow, as well as palm oil, olein, and stearin.^{1,2} 1,2-Dipalmitoyl-3-oleoyl-*rac*-glycerol has been used as a substrate to determine the regioselectivity and substrate specificity of lipases from *Geotrichum* species.³

References

1. Ngo-Duy, C.-C., Destailats, F., Keskitalo, M., *et al.* Triacylglycerols of *Apiaceae* seed oils: Composition and regiodistribution of fatty acids. *Eur. J. Lipid Sci. Technol.* **111**(2), 164-169 (2009).
2. Kallio, H., Yli-Jokipii, K., Kurvinen, J.-P., *et al.* Regioisomerism of triacylglycerols in lard, tallow, yolk, chicken skin, palm oil, palm olein, palm stearin, and a transesterified blend of palm stearin and coconut oil analyzed by tandem mass spectrometry. *J. Agric. Food Chem.* **49**(7), 3363-3369 (2001).
3. Stránský, K., Zarevúcka, M., Kejík, Z., *et al.* Substrate specificity, regioselectivity and hydrolytic activity of lipases activated from *Geotrichum* sp. *Biochem. Eng. J.* **34**(3), 209-216 (2007).

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