

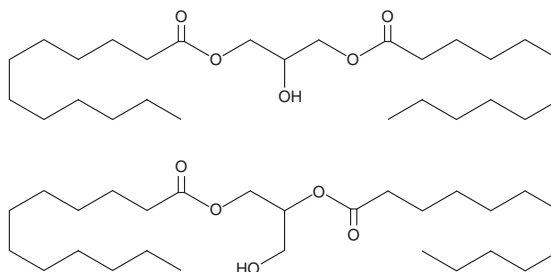
PRODUCT INFORMATION



Dilauroyl-*rac*-glycerol

Item No. 16536

CAS Registry No.: 27638-00-2
Formal Name: diester with 1,2,3-propanetriol dodecanoic acid
Synonyms: Didodecanoylglycerol, Dilaurin
MF: C₂₇H₅₂O₅
FW: 456.7
Purity: ≥98%
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

Dilauroyl-*rac*-glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the dilauroyl-*rac*-glycerol in the solvent of choice. Dilauroyl-*rac*-glycerol is soluble in organic solvents such as ethanol and dimethyl formamide, which should be purged with an inert gas. The solubility of dilauroyl-*rac*-glycerol in these solvents is approximately 1 and 20 mg/ml, respectively.

Description

Dilauroyl-*rac*-glycerol is comprised of a mixture of C12:0 diglyceride racemates with acyl groups attached at either α , α' or α , β -positions. It has been used for the investigation of biological and physicochemical relationships of membrane glycerides.^{1,2} It has also been employed as an industrial surfactant and a building block for synthesis of lipid derivatives.³

References

1. Oo, K.C. and Stumpf, P.K. Some enzymic activities in the germinating oil palm (*Elaeis guineensis*) seedling. *Plant Physiol.* **73**(4), 1028-1032 (1983).
2. Lin, C.Y., Abraham, S., and Smith, S. Acyl specificity in triglyceride synthesis by lactating rat mammary gland. *J. Lipid Res.* **17**(6), 647-656 (1976).
3. Zhong, N., Gui, Z., Huang, J., *et al.* Solvent-free enzymatic synthesis of 1,3-diacylglycerols by direct esterification of glycerol with saturated fatty acids. *Lipids Health Dis.* **12**(1), 1-7 (2013).

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, 10/26/2022

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