

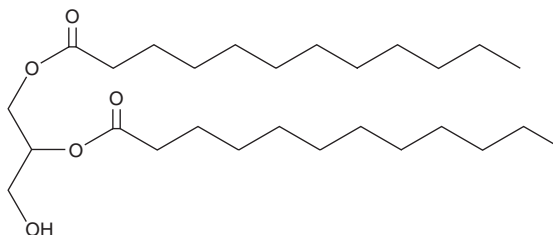
PRODUCT INFORMATION



1,2-Dilauroyl-*rac*-glycerol

Item No. 26846

CAS Registry No.: 17598-94-6
Formal Name: dodecanoic acid, 1-(hydroxymethyl)-
1,2-ethanediyl ester
Synonyms: DG(12:0/12:0/0:0),
1,2-Didodecanoyl-*rac*-glycerol,
1,2-Dilaurin
MF: C₂₇H₅₂O₅
FW: 456.7
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-Dilauroyl-*rac*-glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-dilauroyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1,2-Dilauroyl-*rac*-glycerol is soluble in the organic solvent dimethyl formamide (DMF) at a concentration of approximately 20 mg/ml.

1,2-Dilauroyl-*rac*-glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,2-dilauroyl-*rac*-glycerol should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 1,2-Dilauroyl-*rac*-glycerol has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

1,2-Dilauroyl-*rac*-glycerol is a diacylglycerol that contains lauric acid (Item No. 10006626) at the *sn*-1 and *sn*-2 positions. It has been used as an internal standard for the quantification of diacylglycerols in rat desheathed sciatic nerves.¹ Monomolecular films containing 1,2-dilauroyl-*rac*-glycerol have been used as substrates to measure the effects of surface pressure and pancreatic procolipase and colipase on the activity of porcine pancreatic lipase.²

References

1. Zhu, X. and Eichberg, J. 1,2-Diacylglycerol content and its arachidonyl-containing molecular species are reduced in sciatic nerve from streptozotocin-induced diabetic rats. *J. Neurochem.* **55**(3), 1087-1090 (1990).
2. Wieloch, T., Borgström, B., Piérioni, G., *et al.* Porcine pancreatic procolipase and its trypsin-activated form: Lipid binding and lipase activation on monomolecular films. *FEBS Lett.* **128**(2), 217-220 (1981).

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, 11/08/2022

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