

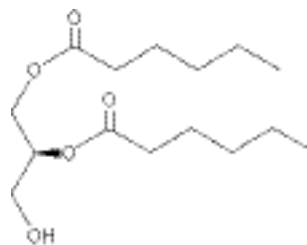
Product Information



1,2-Dihexanoyl-*sn*-glycerol

Catalog No. 10008646

CAS Registry No.: 30403-47-5
Formal Name: 1,2-*bis*(O-hexanoyl)-*sn*-glycerol
MF: C₁₅H₂₈O₅
FW: 288.4
Purity: ≥95%
Stability: ≥1 year at -80°C
Supplied as: A solution in ethanol



Laboratory Procedures

For long term storage, we suggest that 1,2-dihexanoyl-*sn*-glycerol be stored as supplied at -80°C. It will be stable for at least one year.

1,2-Dihexanoyl-*sn*-glycerol is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of 1,2-dihexanoyl-*sn*-glycerol in DMSO is approximately 7 mg/ml and approximately 20 mg/ml in DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 1,2-dihexanoyl-*sn*-glycerol can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 1,2-dihexanoyl-*sn*-glycerol in PBS (pH 7.2) is approximately 250 µg/ml. We do not recommend storing the aqueous solution for more than one day.

1,2-Dihexanoyl-*sn*-glycerol is an analog of the protein kinase C (PKC)-activating second messenger diacylglycerol (DAG). Although the biological activities of 1,2-dihexanoyl-*sn*-glycerol have not been well characterized, it is expected to behave similarly to 1,2-dioctanoyl-*sn*-glycerol (Catalog No. 62225).^{1,2}

References

- Ostroski, M., Tu-Sekine, B., and Raben, D.M. Analysis of a novel diacylglycerol kinase from Dictyostelium discoideum: DGKA. *Biochemistry* **44**, 10199-10207 (2005).
- Doherty, C.M., Tarchala, S.M., Radwanska, E., *et al.* Characterization of two second messenger pathways and their interactions in eliciting the human sperm acrosome reaction. *J. Androl.* **16**, 36-46 (1995).

Related Products

1,2-Didecanoyl-*sn*-glycerol - Cat. No. 62210 • 1,2-Dioctanoyl-*sn*-glycerol - Cat. No. 62225 • 1,2-Dioleoyl-*sn*-glycerol - Cat. No. 62230 • 1-Oleoyl-2-acetyl-*sn*-glycerol - Cat. No. 62600

WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 09/21/2007

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com