

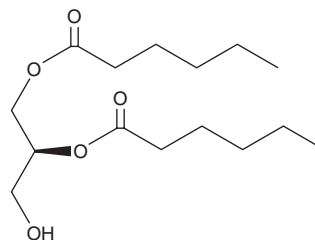
PRODUCT INFORMATION



1,2-Dihexanoyl-*sn*-glycerol

Item No. 10008646

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



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

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 purged with an inert gas can be used. The solubility of 1,2-dihexanoyl-*sn*-glycerol in these solvents is approximately 7 and 20 mg/ml, respectively.

1,2-Dihexanoyl-*sn*-glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 1,2-dihexanoyl-*sn*-glycerol should be diluted with the aqueous buffer of choice. The solubility of 1,2-dihexanoyl-*sn*-glycerol in PBS (pH 7.2) is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

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 (Item No. 62225).^{1,2}

References

1. Ostroski, M., Tu-Sekine, B., and Raben, D.M. Analysis of a novel diacylglycerol kinase from *Dictyostelium discoideum*: DGKA. *Biochemistry* **44(30)**, 10199-10207 (2005).
2. 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(1)**, 36-46 (1995).

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, 03/06/2024

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