

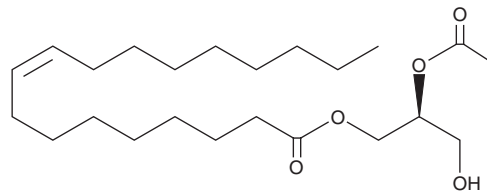
PRODUCT INFORMATION



1-Oleoyl-2-acetyl-sn-glycerol

Item No. 62600

CAS Registry No.: 86390-77-4
Formal Name: 1-O-9Z-octadecenoyl-2-O-acetyl-sn-glycerol
Synonym: OAG
MF: C₂₃H₄₂O₅
FW: 398.6
Purity: ≥95%
Supplied as: A solution in acetonitrile
Storage: -80°C
Stability: ≥1 year



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

Laboratory Procedures

1-Oleoyl-2-acetyl-sn-glycerol (OAG) is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile 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 OAG in these solvents is approximately 20 mg/ml. OAG is also miscible in ethanol.

OAG is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the acetonitrile solution of OAG should be diluted with the aqueous buffer of choice. OAG has a solubility of 1.7 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

OAG is a cell-permeable analog of the PKC-activating second messenger DAG. It activates PKC in platelets, resulting in the phosphorylation of a 40 kDa protein. OAG is metabolized to its corresponding phosphatidic acid (1-oleoyl-2-acetyl-3-phosphoglycerol), most likely through the action of a diglycerol kinase.¹

Reference

1. Nishizuka, Y. The role of protein kinase C in cell surface signal transduction and tumour promotion. *Nature* 308(5961), 693-697 (1984).

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, 02/15/2023

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