

PRODUCT INFORMATION



1-O-hexadecyl-2-Dihomo- γ -Linolenoyl-sn-glycero-3-PC

Item No. 60901

CAS Registry No.: 155575-01-2
Formal Name: 1-O-hexadecyl-2-O-(8Z,11Z,14Z-eicosatrienoyl)-sn-glyceryl-3-phosphorylcholine
Synonyms: 1-O-hexadecyl-2-Dihomo- γ -Linolenoyl-sn-glycero-3-Phosphocholine,
1-O-hexadecyl-2-(8Z,11Z,14Z-eicosatrienoyl)-sn-glycero-Phosphocholine,
1-O-hexadecyl-2-(8Z,11Z,14Z-eicosatrienoyl)-sn-glycero-3-Phosphocholine,
1-O-hexadecyl-2-(8Z,11Z,14Z-eicosatrienoyl)-sn-glyceryl-Phosphocholine,
1-O-hexadecyl-2-O-(8Z,11Z,14Z-eicosatrienoyl)-sn-glyceryl-PC,
PC(O-16:0/20:3)

MF: C₄₄H₈₄NO₇P

FW: 770.1

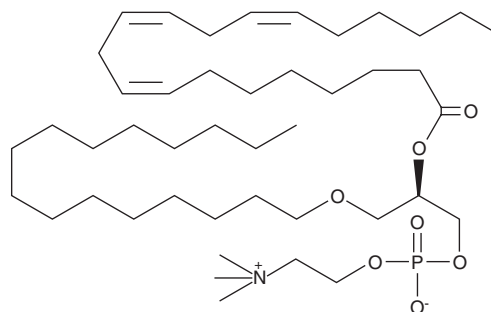
Purity: \geq 98%

Supplied as: A solution in ethanol containing 0.1% BHT

Storage: -20°C

Stability: \geq 1 year

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



Laboratory Procedures

1-O-hexadecyl-2-Dihomo- γ -linolenoyl-sn-glycero-3-PC 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-O-hexadecyl-2-dihomo- γ -linolenoyl-sn-glycero-3-PC in these solvents is approximately 2.5 and 14 mg/ml, respectively.

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. If an organic solvent-free solution of 1-O-hexadecyl-2-dihomo- γ -linolenoyl-sn-glycero-3-PC is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 1-O-hexadecyl-2-dihomo- γ -linolenoyl-sn-glycero-3-PC in PBS (pH 7.2) is approximately 50 μ g/ml. We do not recommend storing the aqueous solution for more than one day.

Description

1-O-hexadecyl-2-Dihomo- γ -linolenoyl-sn-glycero-3-PC is a PAF analog which contains dihomogamma-linolenate at the sn-2 position rather than the acetate moiety found in PAF C-16. The biological activities of 1-O-hexadecyl-2-dihomo- γ -linolenoyl-sn-glycero-3-PC are not known at present, but may be similar to arachidonoyl PAF C-16 (Item No. 60904).

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, 11/15/2021

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