

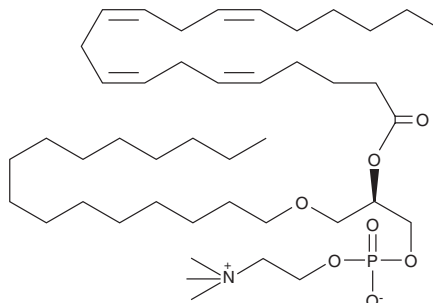
PRODUCT INFORMATION



1-O-hexadecyl-2-Arachidonoyl-*sn*-glycero-3-PC

Item No. 60904

CAS Registry No.: 86288-11-1
Formal Name: 1-O-hexadecyl-2-O-(5Z,8Z,11Z,14Z-eicosatetraenoyl)-*sn*-glyceryl-3-phosphorylcholine
Synonym: Arachidonoyl PAF C-16, 1-O-hexadecyl-2-Arachidonoyl-*sn*-glycero-3-Phosphocholine, PC(O-16:0/20:4)
MF: C₄₄H₈₂NO₇P
FW: 768.2
Purity: ≥95%
Supplied as: A solution in ethanol containing 0.1% BHT
Storage: -20°C
Stability: ≥2 years
Special Conditions: Hygroscopic



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

Laboratory Procedures

1-O-hexadecyl-2-Arachidonoyl-*sn*-glycero-3-PC is supplied as a solution in ethanol containing 0.1% of the antioxidant BHT (2,6-di-*tert*-butyl-4-methyl phenol). 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-arachidonoyl-*sn*-glycero-3-PC in these solvents is approximately 2 and 10 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-arachidonoyl-*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-arachidonoyl-*sn*-glycero-3-PC in PBS (pH 7.2), basic PBS, and acidic PBS is approximately 0.5 mg/ml. Store aqueous solutions of 1-O-hexadecyl-2-arachidonoyl-*sn*-glycero-3-PC on ice and use within 12 hours of preparation. We recommend making a fresh preparation each day.

Description

1-O-hexadecyl-2-Arachidonoyl-*sn*-glycero-3-PC is the product of acylation of lyso-PAF C-16 (Item No. 60906) by a CoA-independent transacylase.^{1,2} It is the most common precursor for formation of PAF C-16 (Item No. 60900) by the remodeling pathway.³

References

1. Kramer, R.M., Patton, G.M., Pritzker, C.R., *et al.* Metabolism of platelet-activating factor in human platelets. Transacylase-mediated synthesis of 1-O-alkyl-2-arachidonoyl-*sn*-glycero-3-phosphocholine. *J. Biol. Chem.* **259(21)**, 13316-13320 (1984).
2. Venable, M.E., Olson, S.C., Nieto, M.L., *et al.* Enzymatic studies of lyso platelet-activating factor acylation in human neutrophils and changes upon stimulation. *J. Biol. Chem.* **268(11)**, 7965-7975 (1993).
3. Prescott, S.M., Zimmerman, G.A., and McIntyre, T.M. Platelet-activating factor. *J. Biol. Chem.* **265(29)**, 17381-17384 (1990).

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, 03/11/2024

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