

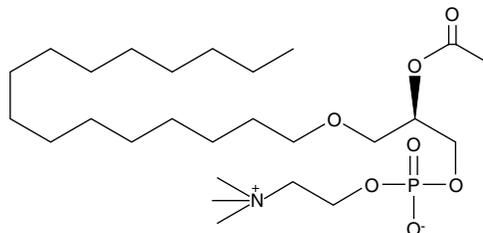
Product Information



PAF C-16

Item No. 60900

CAS Registry No.: 74389-68-7
Formal Name: 1-O-hexadecyl-2-O-acetyl-sn-glyceryl-3-phosphorylcholine
Synonym: 1-O-Hexadecyl-2-acetyl-sn-glycero-3-phosphocholine
MF: C₂₆H₅₄NO₇P
FW: 523.7
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A lyophilized powder



Laboratory Procedures

For long term storage, we suggest that PAF C-16 be stored as supplied at -20°C, preferably in a desiccator. It will be stable for at least one year.

PAF C-16 is supplied as a lyophilized powder. A stock solution may be made by dissolving the PAF C-16 in an organic solvent purged with an inert gas. PAF C-16 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PAF C-16 in these solvents is approximately 10 mg/ml.

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 PAF C-16 can be prepared by directly dissolving the lyophilized powder in aqueous buffers. The solubility of PAF C-16 in PBS, pH 7.2, is approximately 25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

PAF C-16 is a phospholipid synthesized by stimulated endothelial cells, neutrophils, basophils, eosinophils, and macrophages.^{1,2} In addition to activation of platelets, PAF C-16 also activates neutrophils, monocytes, and macrophages. Its receptor mediated biological actions are achieved at concentrations as low as 10⁻¹² M.^{3,4}

References

1. Benveniste, J. Platelet-activating factor, a new mediator of anaphylaxis and immune complex deposition from rabbit and human basophils. *Nature* **249**, 581-582 (1974).
2. Whatley, R.E., Zimmerman, G.A., McIntyre, T.M., *et al.* Lipid metabolism and signal transduction in endothelial cells. *Prog. Lipid Res.* **29**, 45-63 (1990).
3. Hanahan, D.J. and Kumar, R. Platelet activating factor; chemical and biochemical characteristics. *Prog. Lipid Res.* **26**, 1-28 (1987).
4. Prescott, S.M., Zimmerman, G.A., and McIntyre, T.M. Platelet-activating factor. *J. Biol. Chem.* **265**, 17381-17384 (1990).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/60900

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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 Safety Data Sheet, which has been sent via email to 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, 02/05/2014

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