

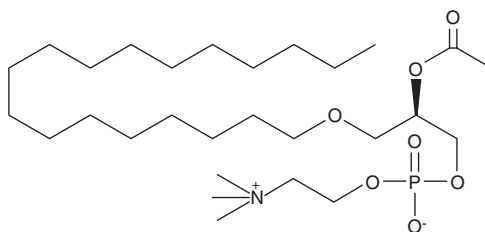
Product Information



PAF C-18

Item No. 60910

CAS Registry No.: 74389-69-8
Formal Name: 1-O-octadecyl-2-O-acetyl-*sn*-glyceryl-3-phosphorylcholine
MF: C₂₈H₅₈NO₇P
FW: 551.7
Purity: ≥97%
Stability: ≥2 years at -20°C
Supplied as: A solution in ethanol



Laboratory Procedures

For long term storage, we suggest that PAF C-18 be stored as supplied at -20°C. It will be stable for at least two years.

PAF C-18 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 PAF C-18 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. If an organic solvent-free solution of PAF C-18 is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of PAF C-18 in PBS (pH 7.2) is approximately 20 mg/ml. We do not recommend storing the aqueous solution for more than one day.

PAF C-18 is a naturally occurring phospholipid produced upon stimulation through two distinct pathways known as the 'remodeling' and 'de novo' pathways.¹ It is less potent than PAF C-16 in the induction of platelet aggregation, but equipotent in activation of guinea pig macrophages.² PAF C-18 induces the release of PGE₂ and TXB₂ from albumin-elicited guinea pig macrophages, and enhances the spreading of plated macrophages.³ Pathological processes involving PAF include necrotizing enterocolitis,⁴ inflammation, asthma, and allergy.⁵

References

1. Prescott, S.M., Zimmerman, G.A., and McIntyre, T.M. Platelet-activating factor. *J. Biol. Chem.* **265**, 17381-17384 (1990).
2. Stewart, A.G. and Grigoriadis, G. Structure-activity relationships for platelet-activating factor (PAF) and analogs reveal differences between PAF receptors on platelets and macrophages. *J. Lipid Mediators* **4**, 299-308 (1991).
3. Hartung, H.-P. Acetyl glyceryl ether phosphorylcholine (platelet-activating factor) mediates heightened metabolic activity in macrophages. Studies on PGE, TXB₂ and O₂ production, spreading, and the influence of calmodulin-inhibitor W-7. *FEBS Lett.* **160**, 209-212 (1983).
4. Wang, H., Tan, X.-D., Qu, X.-W., *et al.* Platelet-activating factor (PAF) up-regulates plasma and tissue PAF-acetylhydrolase activity in the rat: Effect of cycloheximide. *Pediatr. Res.* **42**, 597-603 (1997).
5. Sturk, A., Wouter Ten Cate, J., Hosford, D., *et al.* The synthesis, catabolism, and pathophysiological role of platelet-activating factor. *Adv. Lipid Res.* **23**, 219-276 (1989).

Related Products

PtdIns-(3,4,5)-P₃(1,2-dipalmitoyl)(sodium salt) - Item No. 64920 • Lyso-PAF C-18 -d₄ - Item No. 10010228 • PAF C-18 -d₄ - Item No. 10010229

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

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

MATERIAL 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 Material 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, 11/08/2010