

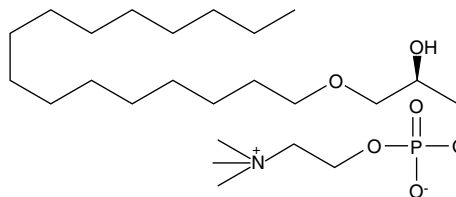
Product Information



Lyso-PAF C-16

Item No. 60906

CAS Registry No.:	52691-62-0
Formal Name:	1-O-hexadecyl- <i>sn</i> -glyceryl-3-phosphorylcholine
Synonyms:	1-O-Hexadecyl- <i>sn</i> -glyceryl-3-phosphocholine
MF:	C ₂₄ H ₅₂ NO ₆ P
FW:	481.7
Purity:	≥98%
Stability:	≥2 years at -20°C
Supplied as:	A lyophilized powder
Misc:	Hygroscopic



Laboratory Procedures

For long term storage, we suggest that Lyso-PAF C-16 be stored as supplied at -20°C. It should be stable for at least two years.

Lyso-PAF C-16 is supplied as a lyophilized powder sealed under vacuum. Prior to performing biological experiments, the lyso-PAF C-16 should be reconstituted in water. The solubility of lyso-PAF C-16 in water is approximately 4.5 mg/ml.

A stock solution may be made by dissolving the lyso-PAF C16 in an organic solvent purged with an inert gas. Lyso-PAF C16 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of lyso-PAF C16 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 lyso-PAF C16 can be prepared by directly dissolving the lyophilized powder in aqueous buffers. The solubility of lyso-PAF C16 in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

PAF C-16 is metabolized to lyso-PAF C16 by the action of PAF acetylhydrolase. Lyso-PAF C16 is recycled in the remodeling pathway of membrane phospholipid biosynthesis.¹⁻³

References

1. Uemura, Y., Lee, T., and Snyder, F. A coenzyme A-independent transacylase is linked to the formation of platelet-activating factor (PAF) by generating the lyso-PAF intermediate in the remodeling pathway. *J. Biol. Chem.* **266**, 8268-8272 (1991).
2. 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).
3. Robinson, M. and Snyder, F. Metabolism of platelet-activating factor by rat alveolar macrophages: lyso-PAF as an obligatory intermediate in the formation of alkylarachidonoyl glycerophosphocholine species. *Biochim. Biophys. Acta* **837**, 52-56 (1985).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/60906

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, 02/14/2014

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