

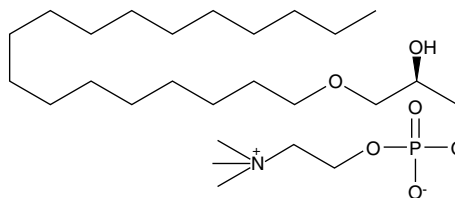
# Product Information



## Lyso-PAF C-18

Item No. 60916

**CAS Registry No:** 74430-89-0  
**Formal Name:** 1-O-octadecyl-*sn*-glyceryl-3-phosphorylcholine  
**MF:** C<sub>26</sub>H<sub>56</sub>NO<sub>6</sub>P  
**FW:** 509.7  
**Purity:** ≥98%  
**Stability:** ≥1 year at -20°C  
**Supplied as:** A lyophilized powder  
**Misc:** Hygroscopic



### Laboratory Procedures

Lyso-PAF C-18 is lysophospholipid which induces PAF biosynthesis. For long term storage, we suggest that lyso-PAF C-18 be stored as supplied at -20°C, preferably in a dessicator. It will be stable for at least one year.

Lyso-PAF C-18 is supplied as a lyophilized powder. Prior to performing biological experiments, the lyophilized powder should be reconstituted in the buffer of choice. Lyso-PAF C-18 is soluble in PBS (pH 7.2) at a concentration of approximately 4.5 mg/ml. The buffer should not contain albumin as the presence of albumin will greatly decrease the amount of PAF produced from lyso-PAF C-18. Sonicate until a clear solution is obtained. Store aqueous solutions of lyso-PAF C-18 on ice and use within 12 hours. We recommend making a fresh preparation each day.

Lyso-PAF C-18 can be formed by either the action of PAF-AH on PAF C-18 or by the action of a CoA-independent transacylase on 1-O-octadecyl-2-acyl-glycerophosphocholine.<sup>1-3</sup> Lyso PAF C-18 is a substrate for either PAF C-18 formation by the remodeling pathway or selective acylation with arachidonic acid by a CoA-independent transacylase.<sup>4,5</sup>

### References

1. Stafforini, D.M., Prescott, S.M., and McIntyre, T.M. Human plasma platelet-activating factor acetylhydrolase. *J. Biol. Chem.* **262**, 4223-4230 (1987).
2. 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).
3. Venable, M.E., Nieto, M.L., Schmitt, J.D., *et al.* Conversion of 1-O-[3H]alkyl-2-arachidonoyl-*sn*-glycero-3-phosphorylcholine to lyso platelet-activating factor by the CoA-independent transacylase in membrane fractions of human neutrophils. *J. Biol. Chem.* **266**, 18691-18698 (1991).
4. Prescott, S.M., Zimmerman, G.A., and McIntyre, T.M. Platelet-activating factor. *J. Biol. Chem.* **265**, 17381-17384 (1990).
5. Venable, M.E., Olson, S.C., Nieto, M.L., *et al.* Enzymatic studies of lyso platelet-activation factor acylation in human neutrophils and changes upon stimulation. *J. Biol. Chem.* **268**, 7965-7975 (1993).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/60916](http://www.caymanchem.com/catalog/60916)

**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 *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, 03/07/2012

### 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](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)