

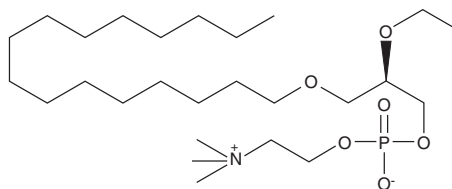
PRODUCT INFORMATION



Lyso-PAF C-16

Item No. 60916

CAS Registry No.: 74430-89-0
Formal Name: 1-O-octadecyl-sn-glycerol-3-phosphorylcholine
MF: $C_{26}H_{56}NO_6P$
FW: 509.7
Purity: $\geq 98\%$
Supplied as: A lyophilized powder
Storage: -20°C
Stability: ≥ 4 years
Special Conditions: Hygroscopic



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

Laboratory Procedures

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.

Description

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.¹⁻³ 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.^{4,5}

References

- Stafforini, D.M., Prescott, S.M., and McIntyre, T.M. Human plasma platelet-activating factor acetylhydrolase. *J. Biol. Chem.* **262**, 4223-4230 (1987).
- 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).
- 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).
- Prescott, S.M., Zimmerman, G.A., and McIntyre, T.M. Platelet-activating factor. *J. Biol. Chem.* **265**, 17381-17384 (1990).
- 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).

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, 05/16/2024

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