

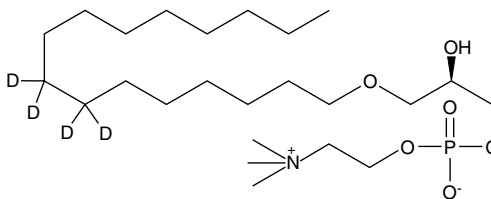
Product Information



Lyso-PAF C-16-d₄

Item No. 360906

CAS Registry No.: 201216-37-7
Formal Name: 1-O-hexadecyl-(7,7,8,8-d₄)-sn-glycero-3-phosphocholine
MF: C₂₄H₅₂D₄NO₆P
FW: 485.7
Chemical Purity: ≥98%
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol



Laboratory Procedures

Lyso-PAF C-16-d₄ contains four deuterium atoms at the 7, 7', 8, and 8' positions of the hexadecyl moiety. It is intended for use as an internal standard for the quantification of lyso-PAF C-16 by GC- or LC-mass spectrometry (MS). For long term storage, we suggest that lyso-PAF C-16-d₄ be stored as supplied at -20°C. It will be stable for at least one year.

Lyso-PAF C-16-d₄ 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 ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of lyso-PAF C-16-d₄ in these solvents is approximately 10 mg/ml.

Lyso-PAF C-16-d₄ is used as an internal standard for the quantification of lyso-PAF C-16-d₄ by stable isotope dilution MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Lyso-PAF C-16 can be formed by either the action of PAF-AH on PAF C-16,¹ or by the action of a CoA-independent transacylase on 1-O-hexadecyl-2-acyl-glycerophosphocholine.^{2,3} Lyso-PAF C-16 is a substrate for either PAF C-16 formation by the remodeling pathway⁴ or selective acylation with arachidonic acid by a CoA-independent transacylase.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).
- Huber, M., Müller, J., Leier, I., *et al.* Metabolism of cysteinyl leukotrienes in monkey and man. *Eur. J. Biochem.* **194**, 309-315 (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).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/360906

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, 09/06/2013

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