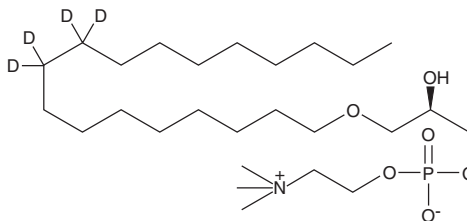


PRODUCT INFORMATION



Lyso-PAF C-18-d₄ Item No. 10010228

Formal Name: 1-O-octadecyl-*sn*-glyceryl-3-phosphorylcholine-9,9,10,10-d₄
MF: C₂₆H₅₂D₄NO₆P
FW: 513.7
Chemical Purity: ≥98% (Lyso-PAF C-18)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Lyso-PAF C-18-d₄ is intended for use as an internal standard for the quantification of lyso-PAF C-18 (Item No. 60916) by GC- or LC-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-18-d₄ is supplied as a crystalline solid. A stock solution may be made by dissolving the lyso-PAF C-18-d₄ in the solvent of choice, which should be purged with an inert gas. Lyso-PAF C-18-d₄ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of lyso-PAF C-18-d₄ in these solvents is approximately 10 mg/ml.

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

1. Stafforini, D.M., Prescott, S.M., and McIntyre, T.M. *J. Biol. Chem.* **262**(9), 4223-4230 (1987).
2. Uemura, Y., Lee, T., and Snyder, F. *J. Biol. Chem.* **266**(13), 8268-8272 (1991).
3. Venable, M.E., Nieto, M.L., Schmitt, J.D., et al. *J. Biol. Chem.* **266**(28), 18691-18698 (1991).
4. Prescott, S.M., Zimmerman, G.A., and McIntyre, T.M. *J. Biol. Chem.* **265**(29), 17381-17384 (1990).
5. Venable, M.E., Olson, S.C., Nieto, M.L., et al. *J. Biol. Chem.* **268**(11), 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, 01/26/2024

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