

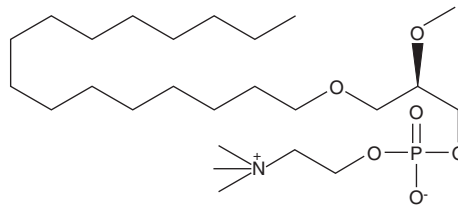
PRODUCT INFORMATION



2-O-methyl PAF C-16

Item No. 60902

CAS Registry No.: 78858-44-3
Formal Name: 1-O-hexadecyl-2-O-methyl-*sn*-glyceryl-3-phosphorylcholine
MF: C₂₅H₅₄NO₆P
FW: 495.7
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

2-O-methyl PAF C-16 is supplied as a solution in ethanol. 2-O-methyl PAF C-16 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 2-O-methyl PAF C-16 should be diluted with the aqueous buffer of choice. The solubility of 2-O-methyl PAF C-16 in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2-O-methyl PAF C-16 is a synthetic PAF analog which contains a methyl group, attached by an ether linkage, at the *sn*-2 position. Although the biological activities of 2-O-methyl PAF C-16 have not been well characterized, the C-18 analog has been shown to mediate a number of biological responses including decreasing plasma membrane fluidity and inhibiting tumor cell invasiveness in embryonic chick heart.¹ In rat astrocytes, the C-18 analog causes the release of large quantities of NO by a pathway involving activation of NOS.²

References

1. van Blitterswijk, W.J., Hilkmann, H., and Storme, G.A. Accumulation of an alkyl lysophospholipid in tumor cell membranes affects membrane fluidity and tumor cell invasion. *Lipids* **22(11)**, 820-823 (1987).
2. Cardile, V., Jiang, X., Renis, M., *et al.* Effects of ether lipid 1-O-octadecyl-2-methoxy-*rac*-glycero-3-phosphocholine and its analogs PAF and CPAF on the release of nitric oxide in primary cultures of rat astrocytes. *Brain Res.* **715(1-2)**, 98-103 (1996).

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, 01/18/2024

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