

PRODUCT INFORMATION

2-O-methyl PAF C-18 (hydrate)

Item No. 60912

Formal Name: 4-hydroxy-7R-methoxy-N,N,N-trimethyl-3,5,9-trioxa-4-phosphaheptacosan-1-aminium, inner salt, 4-oxide, hydrate

Synonyms: Edelfosine, ET-18-OCH₃, 1-O-octadecyl-2-O-methyl-sn-glyceryl-3-phosphorylcholine

MF: C₂₇H₅₈NO₆P • XH₂O

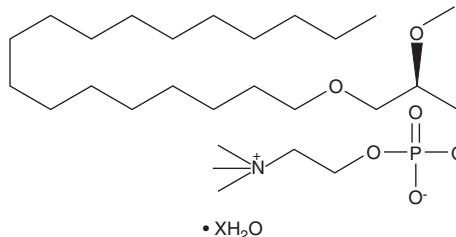
FW: 523.7

Purity: ≥98%

Supplied as: A lyophilized powder

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

2-O-methyl PAF C-18 (hydrate) is supplied as a lyophilized powder. Aqueous solutions of 2-O-methyl PAF C-18 (hydrate) can be prepared by directly dissolving the lyophilized powder in aqueous buffers. The solubility of 2-O-methyl PAF C-18 (hydrate) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2-O-methyl-PAF C-18 is a synthetic PAF analog which contains a methyl group, attached by an ether linkage, at the *sn*-2 position. It 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, 2-O-methyl PAF C-18 causes the release of large quantities of nitric oxide by a pathway involving activation of nitric oxide synthase.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 Research* **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, 11/02/2022

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