

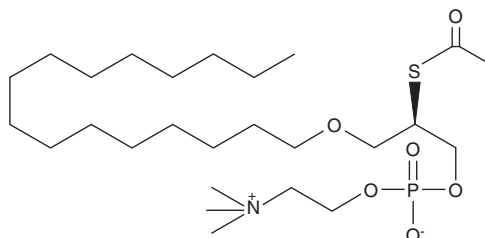
PRODUCT INFORMATION



2-thio PAF

Item No. 60945

CAS Registry No.: 96801-55-7
Formal Name: 1-O-hexadecyl-2-deoxy-2-thio-S-acetyl-sn-glyceryl-3-phosphorylcholine
MF: C₂₆H₅₄NO₆PS
FW: 539.8
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -80°C
Stability: ≥2 years



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

Laboratory Procedures

2-thio PAF 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 2-thio PAF in these solvents is approximately 25 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of 2-thio PAF is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 2-thio PAF in PBS (pH 7.2) is approximately 50 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2-thio PAF is an isosteric analog of PAF. It is also a synthetic analog of PAF which contains an acetyl group attached by a thioester bond at the sn-2 position. It functions as a PAF receptor agonist, and its potency is comparable to PAF C-18 in rabbit platelet aggregation, and to PAF C-16 in the activation of guinea pig macrophages.¹ 2-thio PAF is a substrate for lysophospholipase II and for PAF-AH, and is the substrate used in Cayman's PAF-AH Assay Kit (Item No. 760901).²

References

1. Stewart, A.G. and Grigoriadis, G. Structure-activity relationships for platelet-activating factor (PAF) and analogs reveal differences between PAF receptors on platelets and macrophages. *J. Lipid Mediat.* **4**(3), 299-308 (1991).
2. Aarsman, A.J., Neys, F.W., and Van den Bosch, H. Catabolism of platelet-activating factor and its acyl analog. Differentiation of the activities of lysophospholipase and platelet-activating-factor acetylhydrolase. *Eur. J. Biochem.* **200**(1), 187-193 (1991).

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, 03/11/2024

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