

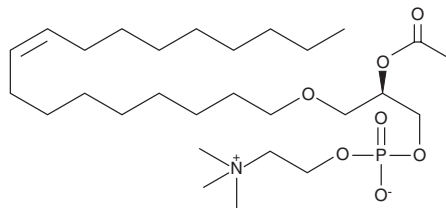
# PRODUCT INFORMATION



## PAF C-18:1

Item No. 18779

**CAS Registry No.:** 85966-90-1  
**Formal Name:** 7R-(acetyloxy)-4-hydroxy-N,N,N-trimethyl-3,5,9-trioxa-4-phosphaheptacos-18Z-en-1-aminium, 4-oxide inner salt  
**Synonyms:** 1-Oleyl-2-acetyl-*sn*-glycero-3-phosphorylcholine, Platelet-activating Factor C-18:1  
**MF:** C<sub>28</sub>H<sub>56</sub>NO<sub>7</sub>P  
**FW:** 549.7  
**Purity:** ≥95%  
**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

PAF C-18:1 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 PAF C-18:1 in these solvents is approximately 10 mg/ml. PAF C-18:1 is also soluble in water at a concentration of 20 mg/ml.

PAF C-18:1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of PAF C-18:1 should be diluted with the aqueous buffer of choice. The solubility of PAF C-18:1 in PBS (pH 7.2) is approximately 20 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

PAF C-18:1 is a naturally occurring phospholipid produced by cells upon stimulation and plays a role in the establishment and maintenance of the inflammatory response. It is less potent than PAF C-16 (Item No. 60900) and PAF C-18 (Item No. 60910) in the induction of neutrophil chemotaxis, but is equipotent to PAF C-16 and PAF C-18 in promoting eosinophil migration.<sup>1,2</sup> PAF C-18:1 activates the PAF receptor and has been used in antibody binding experiments to determine the importance of an acyl linkage at the *sn*-2 position for recognition at this receptor.<sup>3</sup>

### References

1. Carolan, E.J. and Casale, T.B. Degree of platelet activating factor-induced neutrophil migration is dependent upon the molecular species. *J. Immunol.* **145**(8), 2561-2565 (1990).
2. Erger, R.A. and Casale, T.B. Eosinophil migration in response to three molecular species of platelet activating factor. *Inflamm. Res.* **45**(6), 265-267 (1996).
3. Smal, M.A., Baldo, B.A., and Harle, D.G. The specificity of the binding of platelet activating factor (PAF) to anti-PAF antibodies. *J. Mol. Recognit.* **3**(4), 169-173 (1990).

#### 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, 03/07/2025

#### 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](http://WWW.CAYMANCHEM.COM)