

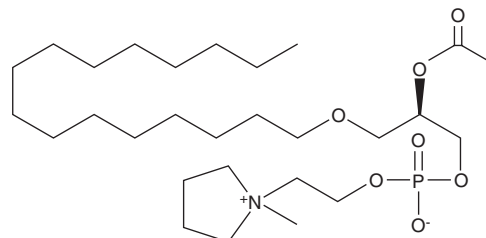
# PRODUCT INFORMATION



## Pyrrolidino PAF C-16

Item No. 60909

**CAS Registry No.:** 91021-63-5  
**Formal Name:** 1-O-hexadecyl-2-O-acetyl-sn-glyceryl-3-phosphoryl-N-methyl-pyrrolidinium ethanol  
**MF:** C<sub>28</sub>H<sub>56</sub>NO<sub>7</sub>P  
**FW:** 549.7  
**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

Pyrrolidino PAF C-16 is supplied as a lyophilized powder. Aqueous solutions of pyrrolidino PAF C-16 can be prepared by directly dissolving the lyophilized powder in aqueous buffers. The solubility of pyrrolidino PAF C-16 in PBS (pH 7.2) is approximately 25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Pyrrolidino PAF C-16 is a synthetic analog of PAF C-16. It is a potent mediator of PAF-like activities, including hypotension<sup>1,2</sup> and platelet aggregation, where it is approximately 3-10 times more potent than PAF itself.<sup>1,2</sup>

### References

1. Shigenobu, K., Masuda, Y., Tanaka, Y., *et al.* Platelet activating factor analogues: Lack of correlation between their activities to produce hypotension and endothelium-mediated vasodilation. *J. Pharmacobiodyn.* **8(2)**, 128-133 (1985).
2. Kasuya, Y., Masuda, Y., and Shigenobu, K. Hypotensive actions of some analogues of platelet-activating factor (PAF) with higher potencies than natural PAF. *Can. J. Physiol. Pharmacol.* **62(4)**, 457-459 (1984).

#### 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, 11/17/2025

#### 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