

# PRODUCT INFORMATION



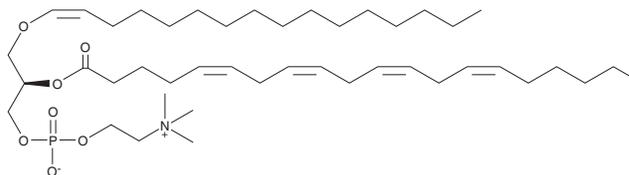
## 1-1(Z)-Hexadecenyl-2-Arachidonoyl-*sn*-glycero-3-PC

Item No. 34182

**CAS Registry No.:** 84460-45-7  
**Formal Name:** (7R)-7-[[[(1Z)-1-hexadecen-1-yloxy]methyl]-4-hydroxy-N,N,N-trimethyl-9-oxo-3,5,8-trioxa-4-phosphaoctacos-13Z,16Z,19Z,22Z-tetraen-1-aminium, inner salt, 4-oxide

**Synonyms:** 1-1(Z)-Hexadecenyl-2-Arachidonoyl-*sn*-glycero-3-Phosphocholine,  
1-1(Z)-Hexadecenyl-2-Arachidonoyl-*sn*-glycero-3-Phosphatidylcholine,  
16:0p/20:4-PC, PC(P-16:0/20:4),  
C16(plasm)-20:4-PC, PlgPC 16:0/20:4

**MF:** C<sub>44</sub>H<sub>80</sub>NO<sub>7</sub>P  
**FW:** 766.1  
**Purity:** ≥95%  
**Supplied as:** A 1 mg/ml solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥1 year



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

### Description

1-1(Z)-Hexadecenyl-2-arachidonoyl-*sn*-glycero-3-PC is a plasmalogen that contains 1(Z)-hexadecenoic acid and arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) at the *sn*-1 and *sn*-2 positions, respectively. Serum levels of 1-1(Z)-hexadecenyl-2-arachidonoyl-*sn*-glycero-3-PC are decreased in dogs with myxomatous mitral valve disease when fed a diet including medium chain triglycerides, fish oil, carnitine precursors, magnesium, and antioxidants.<sup>1</sup>

### Reference

1. Li, Q., Laflamme, D.P., and Bauer, J.E. Serum untargeted metabolomic changes in response to diet intervention in dogs with preclinical myxomatous mitral valve disease. *PLoS One* **15(6)**, e0234404 (2020).

**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, 01/05/2026

### 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