

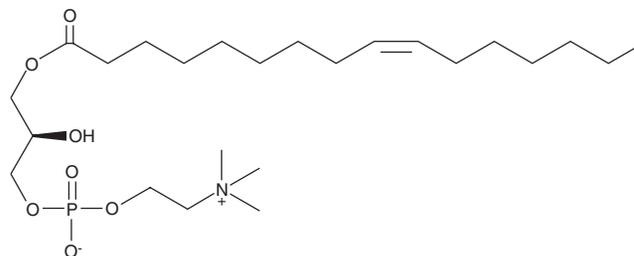
# PRODUCT INFORMATION



## 1-Palmitoleoyl-2-hydroxy-*sn*-glycero-3-PC

Item No. 40522

**CAS Registry No.:** 76790-27-7  
**Formal Name:** (7R,18Z)-4,7-dihydroxy-N,N,N-trimethyl-10-oxo-3,5,9-trioxa-4-phosphapentacos-18-en-1-aminium, 4-oxide, inner salt  
**Synonyms:** 1-Hexadecenoyl-2-hydroxy-*sn*-glycero-3-PC, LPC(16:1), 16:1 Lyso PC, PC(16:1/0:0)  
**MF:** C<sub>24</sub>H<sub>48</sub>NO<sub>7</sub>P  
**FW:** 493.6  
**Purity:** ≥98% (as a 9:1 mixture of the 1-acyl and 2-acyl)  
**Supplied as:** A 10 mg/ml 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.

### Description

1-Palmitoleoyl-2-hydroxy-*sn*-glycero-3-PC is a lysophospholipid. It increases cAMP production in human pancreatic  $\beta$  cells, an effect that can be blocked by the GPR55 antagonist CID-16020046 (Item No. 15247) or the GPR119 antagonist C8.<sup>1</sup> 1-Palmitoleoyl-2-hydroxy-*sn*-glycero-3-PC increases glucose-stimulated insulin secretion in the same cell type, an effect that can be reduced by CID-16020046 or C8.

### Reference

1. Szustak, M., Korkus, E., Madaj, R., *et al.* Lysophosphatidylcholines enriched with cis and trans palmitoleic acid regulate insulin secretion via GPR119 receptor. *ACS Med. Chem. Lett.* **15(2)**, 197-204 (2024).

#### 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, 11/18/2025

#### 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