

# PRODUCT INFORMATION



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

Item No. 33331

**CAS Registry No.:** 50930-23-9  
**Formal Name:** 4,7-dihydroxy-N,N,N-trimethyl-10-oxo-3,5,9-trioxa-4-phosphahehexosan-1-aminium, 4-oxide, inner salt

**Synonyms:** LPC(17:0), 17:0 Lyso-PC, PC(17:0/0:0), 17:0/0:0-PC

**MF:** C<sub>25</sub>H<sub>52</sub>NO<sub>7</sub>P

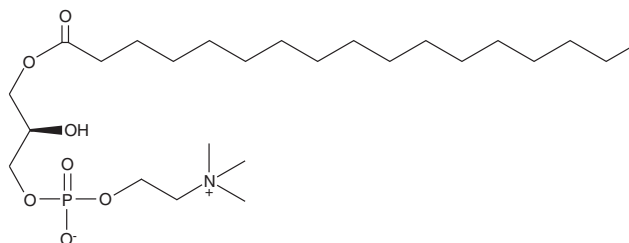
**FW:** 509.7

**Purity:** ≥95%

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥4 years



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

### Laboratory Procedures

1-Heptadecanoyl-2-hydroxy-*sn*-glycero-3-PC is supplied as a crystalline solid. A stock solution may be made by dissolving the 1-heptadecanoyl-2-hydroxy-*sn*-glycero-3-PC in the solvent of choice, which should be purged with an inert gas. 1-Heptadecanoyl-2-hydroxy-*sn*-glycero-3-PC is soluble in the organic solvent ethanol at a concentration of approximately 25 mg/ml.

### Description

1-Heptadecanoyl-2-hydroxy-*sn*-glycero-3-PC is a lysophospholipid that contains the odd-chain saturated fatty acid heptadecanoic acid (Item No. 19722) at the *sn*-1 position. Plasma levels of 1-heptadecanoyl-2-hydroxy-*sn*-glycero-3-PC are increased in a mouse model of Alzheimer's disease induced by intracerebroventricular administration of amyloid- $\beta$  1-42 (A $\beta$ 42; Item No. 20574).<sup>1</sup>

### Reference

1. Gong, Y., Liu, Y., Zhou, L., *et al.* A UHPLC-TOF/MS method based metabonomic study of total ginsenosides effects on Alzheimer disease mouse model. *J. Pharm. Biomed. Anal.* **115**, 174-182 (2015).

#### 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, 04/15/2024

#### 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