

PRODUCT INFORMATION



1-Stearoyl-2-hydroxy-*sn*-glycero-3-PC

Item No. 15107

CAS Registry No.: 19420-57-6
Formal Name: 1-stearoyl-2-lyso-*sn*-glycero-3-phosphatidylcholine
Synonyms: C18:0-PC, Lyso-PC,
1-Octadecanoyl-2-hydroxy-*sn*-glycero-3-phosphatidylcholine,
PC(18:0/0:0), 1-Stearoyl-2-hydroxy-*sn*-glycero-3-Phosphocholine,
1-Stearoyl-2-lyso-*sn*-glycero-3-PC

MF: C₂₆H₅₄NO₇P

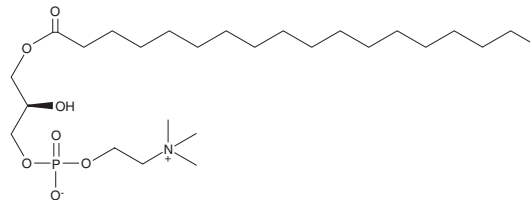
FW: 523.7

Purity: ≥98%

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

Description

1-Stearoyl-2-hydroxy-*sn*-glycero-3-PC is a saturated lysophosphatidylcholine. It increases IL-1 β production in isolated human blood monocytes when used at a concentration of 12.5 μ M.¹ Plasma levels of 1-stearoyl-2-hydroxy-*sn*-glycero-3-PC are decreased in patients with lung cancer.²

References

1. Liu-Wu, Y., Hurt-Camejo, E., and Wiklund, O. Lysophosphatidylcholine induces the production of IL-1 β by human monocytes. *Atherosclerosis* **137**(2), 351-357 (1998).
2. Dong, J., Xiaoming, C., Zhao, L., *et al.* Lysophosphatidylcholine profiling of plasma: Discrimination of isomers and discovery of lung cancer biomarkers. *Metabolomics* **6**(4), 478-488 (2010).

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, 11/15/2022

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