

PRODUCT INFORMATION



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

Item No. 17153

CAS Registry No.: 22252-07-9

Formal Name: (7R,18Z,21Z)-4,7-dihydroxy-N,N,N-trimethyl-10-oxo-3,5,9-trioxa-4-phosphaheptacos-18,21-dien-1-aminium, 4-oxide, inner salt

Synonyms: LGPC, 1-Linoleoyl-2-hydroxy-*sn*-glycero-3-Phosphocholine, 1-Linoleoyl-2-hydroxy-*sn*-glycero-3-Phosphatidylcholine, PC(18:2/0:0), 18:2/0:0-PC

MF: C₂₆H₅₀NO₇P

FW: 519.7

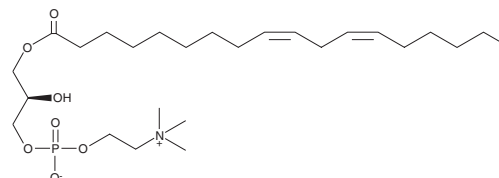
Purity: ≥95%

Supplied as: A 5 mg/ml solution in chloroform

Storage: -20°C

Stability: ≥2 years

Special Conditions: Hygroscopic



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

Laboratory Procedures

1-Linoleoyl-2-hydroxy-*sn*-glycero-3-PC (LGPC) is supplied as a solution in chloroform. LGPC is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the chloroform solution of LGPC should be diluted with the aqueous buffer of choice. The solubility of LGPC in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

LGPC is a lysophospholipid containing linoleic acid (Item Nos. 90150 | 90150.1 | 21909) at the *sn*-1 position that has been found in mouse heart, lung, liver, spleen, kidney, plasma, and serum.¹ Serum levels of LGPC decrease with increasing insulin resistance and dysglycemia in humans.²

References

- Okudaira, M., Inoue, A., Shuto, A., *et al.* Separation and quantification of 2-acyl-1-lysophospholipids and 1-acyl-2-lysophospholipids in biological samples by LC-MS/MS. *J. Lipid Res.* **55**(10), 2178-2192 (2014).
- Gall, W.E., Beebe, K., Lawton, K.A., *et al.* α-Hydroxybutyrate is an early biomarker of insulin resistance and glucose intolerance in a nondiabetic population. *PLoS One* **5**(5), e10883 (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, 11/13/2025

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