

PRODUCT INFORMATION

1-Myristoyl-d₇-2-hydroxy-*sn*-glycero-3-PC

Item No. 9004418

Formal Name: (R)-2-hydroxy-3-((tetradecanoyl-12,12,13,13,14,14,14-d₇)oxy)propyl (2-(trimethylammonio)ethyl) phosphate

Synonyms: 14:0-d₇ Lyso-PC, 1-Myristoyl-d₇-*sn*-glycero-3-Phosphatidylcholine

MF: C₂₂H₃₉D₇NO₇P

FW: 474.6

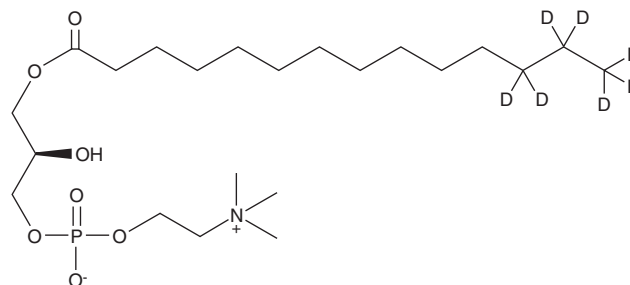
Chemical Purity: ≥95% (1-Myristoyl-2-hydroxy-*sn*-glycero-3-PC)

Deuterium Incorporation: ≥99% deuterated forms (d₁-d₇); ≤1% d₀

Supplied as: A 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-Myristoyl-d₇-2-hydroxy-*sn*-glycero-3-PC is intended for use as an internal standard for the quantification of 1-myristoyl-2-hydroxy-*sn*-glycero-3-PC (Item No. 20958) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Description

1-Myristoyl-d₇-2-hydroxy-*sn*-glycero-3-PC is intended for use as an internal standard for the quantification of 1-myristoyl-2-hydroxy-*sn*-glycero-3-PC (Item No. 20958) by GC- or LC-MS. 1-Myristoyl-2-hydroxy-*sn*-glycero-3-PC is a saturated 14:0 lysophosphatidylcholine that consists of one chain of myristic acid (Item No. 13351) at the *sn*-1 position. It is formed in plasma by lecithin:cholesterol acyltransferase (LCAT) and has a role in lipid signaling by acting on lysophospholipid receptors.¹

Reference

1. Soupene, E., Fyrst, H., and Kuypers, F.A. Mammalian acyl-CoA:lysophosphatidylcholine acyltransferase enzymes. *Proc. Natl. Acad. Sci. USA* **105**(1), 88-93 (2008).

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, 02/03/2025

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