

PRODUCT INFORMATION



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

Item No. 35675

Formal Name: 4-hydroxy-N,N,N-trimethyl-7R-[[[(1-oxotetradecyl-12,12,13,13,14,14,14-d₇)oxy]methyl]-4-oxide, 3,5,8-trioxa-4-phosphahexacos-17Z,20Z-dien-1-aminium, inner salt

Synonyms: 1-Myristoyl-d₇-2-Linoleoyl-*sn*-glycero-3-Phosphatidylcholine, 1-Myristoyl-d₇-2-Linoleoyl-*sn*-glycero-3-Phosphocholine, PC(14:0-d₇/18:2), 14:0-d₇/18:2-PC

MF: C₄₀H₆₉D₇NO₈P

FW: 737.1

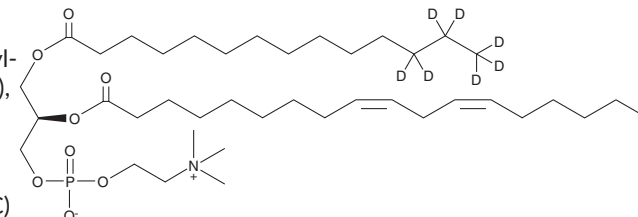
Chemical Purity: ≥98% (1-Myristoyl-2-Linoleoyl-*sn*-glycero-3-PC)

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

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥1 year



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

Laboratory Procedures

1-Myristoyl-d₇-2-linoleoyl-*sn*-glycero-3-PC is intended for use as an internal standard for the quantification of 1-myristoyl-2-linoleoyl-*sn*-glycero-3-PC 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-2-linoleoyl-*sn*-glycero-3-PC is a phospholipid that contains myristic acid (Item No. 13351) and linoleic acid (Item Nos. 90150 | 90150.1 | 21909) at the *sn*-1 and *sn*-2 positions, respectively. It has been found in human plasma.¹

Reference

1. Uhl, O., Glaser, C.B., Demmelmair, H., *et al.* Reversed phase LC/MS/MS method for targeted quantification of glycerophospholipid molecular species in plasma. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **879(30)**, 3556-3564 (2011).

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, 02/22/2024

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