

# PRODUCT INFORMATION



## 1-Palmitoyl-2-13(S)-HODE-*sn*-glycero-3-PC

Item No. 41445

CAS Registry No.: 161923-56-4

Formal Name: (7R,17Z,19E,21S)-4,21-dihydroxy-N,N,N-trimethyl-9-oxo-7-[[[(1-oxohexadecyl)oxy]methyl]-3,5,8-trioxa-4-phosphahexacos-17,19-dien-1-aminium, inner salt, 4-oxide

Synonyms: 1-Hexadecanoyl-2-13(S)-HODE-*sn*-glycero-3-Phosphocholine, 13(S)-Hydroxyoctadecadienoic Acid-PLPC, 1-Palmitoyl-2-13(S)-HODE-*sn*-glycero-3-Phosphocholine, 1-Palmitoyl-2-(13-hydroxy-*cis*-9,*trans*-11-octadecadienoyl)-L-3-phosphatidylcholine

MF:  $C_{42}H_{80}NO_9P$

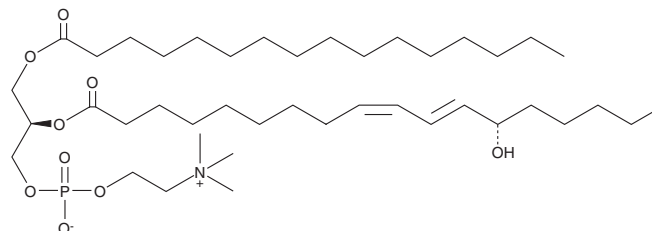
FW: 774.1

Purity:  $\geq 95\%$

Supplied as: A solution in ethanol

Storage:  $-80^{\circ}\text{C}$

Stability:  $\geq 1$  year



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

### Description

1-Palmitoyl-2-13(S)-HODE-*sn*-glycero-3-PC is an oxidized phospholipid that contains palmitic acid (Item No. 10006627) and 13(S)-HODE (Item No. 38610) at the *sn*-1 and *sn*-2 positions, respectively. It has been used to detect the activity of glutathione peroxidase 4 (GPX4), also known as phospholipid hydroperoxide glutathione peroxidase (PhGPx), using 1-palmitoyl-2-13(S)-HpODE-*sn*-glycero-3-PC (Item No. 31976) as the substrate.<sup>1</sup>

### Reference

1. Bao, Y., Chambers, S.J., and Williamson, G. Direct separation of hydroperoxy- and hydroxy-phosphatidylcholine derivatives: Application to the assay of phospholipid hydroperoxide glutathione peroxidase. *Anal. Biochem.* **224**(1), 395-399 (1995).

#### 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, 09/09/2024

#### 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