

PRODUCT INFORMATION



1-Palmitoyl-2-(±)17(18)-EpETE-*sn*-glycero-3-PC

Item No. 24383

Formal Name: (R)-2-(((5Z,8Z,11Z,14Z)-16-((2R,3S)-3-ethyloxiran-2-yl)hexadeca-5,8,11,14-tetraenoyloxy)-3-(palmitoyloxy)propyl (2-(trimethylammonio)ethyl) phosphate

Synonyms: 1-Palmitoyl-2-(17,18-EpETE)-PC, 1-Palmitoyl-2-(17,18-EpETE)-*sn*-glycero-3-PC, 1-Palmitoyl-2-(±)17(18)-EpETE-*sn*-glycero-3-Phosphatidylcholine, 16:0/17,18-EpETE-PC

MF: C₄₄H₇₈NO₉P

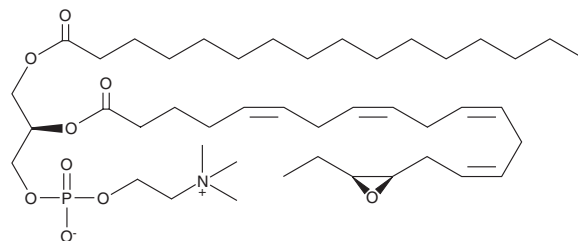
FW: 796.1

Purity: ≥95%

Supplied as: A 100 µg/ml solution in ethanol

Storage: -20°C

Stability: ≥2 years



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

Description

1-Palmitoyl-2-(±)17(18)-EpETE-*sn*-glycero-3-PC is an ω-3 epoxidated eicosapentaenoic acid- and palmitic acid-containing form of phosphatidylcholine.¹ 1-Palmitoyl-2-(±)17(18)-EpETE-*sn*-glycero-3-PC is hydrolyzed by type II platelet activating factor acetylhydrolase (PAF-AH2), but not mutant PAF-AH2 (S234C), to release (±)17(18)-EpETE (Item No. 50861).

Reference

1. Shimanaka, Y., Kono, N., Taketomi, Y., *et al.* Omega-3 fatty acid epoxides are autocrine mediators that control the magnitude of IgE-mediated mast cell activation. *Nat. Med.* **23(11)**, 1287-1297 (2017).

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, 01/29/2026

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