

PRODUCT INFORMATION

1-Palmitoyl-2-linoleoyl-*sn*-glycero-3-phosphoethanolamine Epoxides MaxSpec® LC-MS Mixture

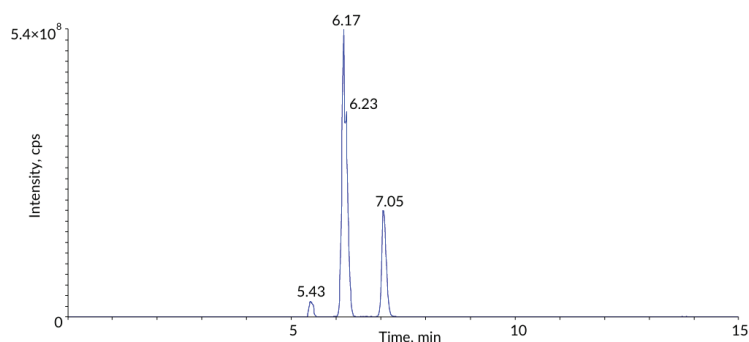
Item No. 45891

Supplied as: A 100 µg/ml mixture in ethanol; in a deactivated glass ampule
Fill Volume: >1 ml
Storage: -20°C (Store upright and unopened)
Stability: ≥2 years
Special Conditions: Warm to room temperature prior to opening. Light sensitive.

Description

1-Palmitoyl-2-linoleoyl-*sn*-glycero-3-phosphoethanolamine Epoxides MaxSpec® LC-MS Mixture (Item No. 45891) is a mixture containing epoxides resulting from mCPBA-mediated epoxidation. It also contains the unoxidized phospholipid, 1-palmitoyl-2-linoleoyl-*sn*-glycero-3-phosphoethanolamine (1,2-PLPE; Item No. 10007072). It may be used as a reference standard for LC-MS to identify phospholipids containing esterified oxylipins in samples derived from a natural source. The mixture is provided as a 1 ml ethanolic solution at 100 µg/ml in an amber ampule in which the headspace has been purged with argon to prevent further lipid oxidation. After opening, it is recommended that the mixture be transferred immediately to a glass screw cap vial to prevent solvent evaporation and stored at -20°C, avoiding multiple freeze/thaw cycles. Ongoing stability testing is performed throughout the shelf life of the product. *NOTE: The amount of solution added to the vial is in excess of the listed amount. Follow recommended storage and handling conditions to maintain product quality.*

Contents



Item No.	Item Name	Retention Time (min)
45889	1-Palmitoyl-2-(9,10;12,13-diepoxyoctadecanoyl)- <i>sn</i> -glycero-3-phosphoethanolamine	5.43
N/A	1-Palmitoyl-2-(9,10-epoxy-12-octadecenoyl)- <i>sn</i> -glycero-3-phosphoethanolamine	6.23
N/A	1-Palmitoyl-2-(12,13-epoxy-9-octadecenoyl)- <i>sn</i> -glycero-3-phosphoethanolamine	6.17
10007072	1-Palmitoyl-2-linoleoyl- <i>sn</i> -glycero-3-phosphoethanolamine	7.05

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, 06/02/2026

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