

PRODUCT INFORMATION

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

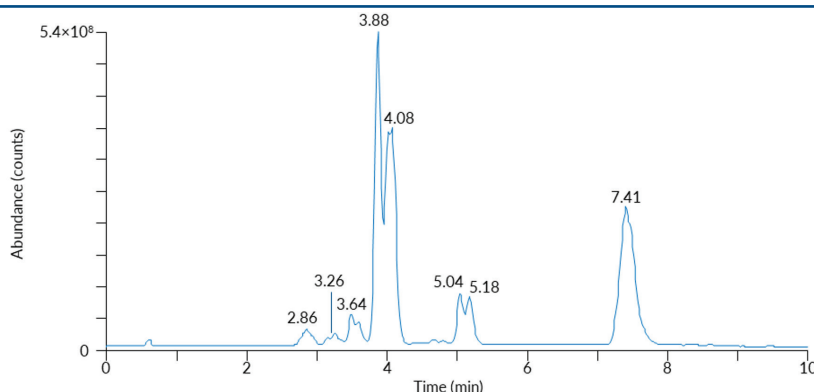
Item No. 46039

Synonym: 1,2-PLPE Vicinal Diol
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 Vicinal Diols MaxSpec® LC-MS Mixture (Item No. 46039) is a mixture containing vicinal diols and epoxides of the alkenes present in 1-palmitoyl-2-linoleoyl-*sn*-glycero-3-phosphoethanolamine (Item No. 10007072). The mixture also contains the unoxidized phospholipid precursor. 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)
N/A	1-Palmitoyl-2-(9,10,12,13-tetrahydroxyoctadecanoyl)- <i>sn</i> -glycero-3-phosphoethanolamine	2.86
N/A	1-Palmitoyl-2-(9,10-epoxy-12,13-dihydroxyoctadecanoyl)- <i>sn</i> -glycero-3-phosphoethanolamine	3.26-3.64
N/A	1-Palmitoyl-2-(12,13-epoxy-9,10-dihydroxyoctadecanoyl)- <i>sn</i> -glycero-3-phosphoethanolamine	3.26-3.64
N/A	1-Palmitoyl-2-(12,13-dihydroxy-9-octadecenoyl)- <i>sn</i> -glycero-3-phosphoethanolamine	3.88
N/A	1-Palmitoyl-2-(9,10-dihydroxy-12-octadecenoyl)- <i>sn</i> -glycero-3-phosphoethanolamine	4.08
45889	1-Palmitoyl-2-(9,10,12,13-diepoxyoctadecanoyl)- <i>sn</i> -glycero-3-phosphoethanolamine	4.02
N/A	1-Palmitoyl-2-(9,10-epoxy-12-octadecenoyl)- <i>sn</i> -glycero-3-phosphoethanolamine	5.04-5.18
N/A	1-Palmitoyl-2-(12,13-epoxy-9-octadecenoyl)- <i>sn</i> -glycero-3-phosphoethanolamine	5.04-5.18
10007072	1-Palmitoyl-2-linoleoyl- <i>sn</i> -glycero-3-phosphoethanolamine	7.41

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, 07/13/2026

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