

PRODUCT INFORMATION

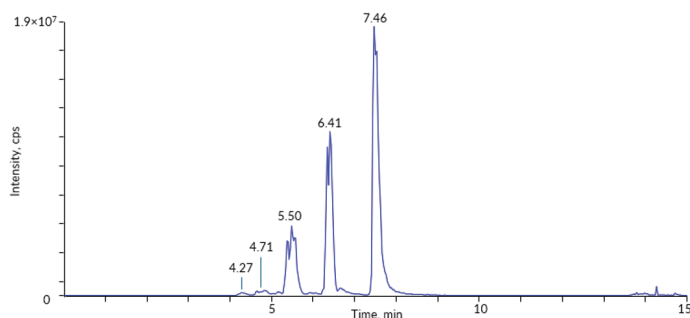
1-Palmitoyl-2-linoleoyl-*sn*-glycero-3-phosphocholine Vicinal Diols MaxSpec® LC-MS Mixture Item No. 45890

Synonym: 1,2-PLPC Vicinal Diols
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-phosphocholine Vicinal Diols MaxSpec® LC-MS Mixture (Item No. 45890) is a mixture containing vicinal diols resulting from oxidation of the alkenes present in 1-palmitoyl-2-linoleoyl-*sn*-glycero-3-phosphocholine (1,2-PLPC; Item No. 20962). 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-phosphocholine	4.27
N/A	1-Palmitoyl-2-(9,10-epoxy-12,13-dihydroxyoctadecanoyl)- <i>sn</i> -glycero-3-phosphocholine	4.66
N/A	1-Palmitoyl-2-(12,13-epoxy-9,10-dihydroxyoctadecanoyl)- <i>sn</i> -glycero-3-phosphocholine	4.83
N/A	1-Palmitoyl-2-(12,13-dihydroxy-9-octadecenoyl)- <i>sn</i> -glycero-3-phosphocholine	5.38
N/A	1-Palmitoyl-2-(9,10-dihydroxy-12-octadecenoyl)- <i>sn</i> -glycero-3-phosphocholine	5.55
45780	1-Palmitoyl-2-(9,10;12,13-diepoxyoctadecanoyl)- <i>sn</i> -glycero-3-phosphocholine	5.55
N/A	1-Palmitoyl-2-(9,10-epoxy-12-octadecenoyl)- <i>sn</i> -glycero-3-phosphocholine	6.32
N/A	1-Palmitoyl-2-(12,13-epoxy-9-octadecenoyl)- <i>sn</i> -glycero-3-phosphocholine	6.41
20962	1-Palmitoyl-2-linoleoyl- <i>sn</i> -glycero-3-phosphocholine	7.46

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, 06/24/2026

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