

PRODUCT INFORMATION



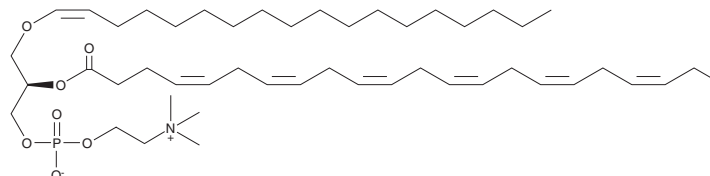
1-1(Z)-Octadecenyl-2-Docosahexaenoyl-*sn*-glycero-3-PC

Item No. 37134

CAS Registry No.: 799268-64-7

Formal Name: 4-hydroxy-N,N,N-trimethyl-7R-[[[(1Z)-1-octadecen-1-yloxy]methyl]-9-oxo-3,5,8-trioxa-4-phosphatriaconta-12Z,15Z,18Z,21Z,24Z,27Z-hexaen-1-

Synonyms: C18(plasm)-22:6-PC,
1-1(Z)-Octadecenyl-2-Docosahexaenoyl-*sn*-glycero-3-Phosphatidylcholine,
1-1(Z)-Octadecenyl-2-Docosahexaenoyl-*sn*-glycero-3-Phosphocholine,
PC(P-18:0/22:6), 18:0p/22:6-PC



MF: C₄₈H₈₄NO₇P

FW: 818.2

Purity: ≥95%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥1 year

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

Description

1-1(Z)-Octadecenyl-2-docosahexaenoyl-*sn*-glycero-3-PC is a plasmalogen that contains 1(Z)-octadecenoic acid and docosahexaenoic acid (DHA; Item Nos. 90310 | 17950) at the *sn*-1 and *sn*-2 positions, respectively. It has been found in bovine retina and fish viscera.^{1,2}

References

- Berdeaux, O., Juaneda, P., Martine, L., *et al.* Identification and quantification of phosphatidylcholines containing very-long-chain polyunsaturated fatty acid in bovine and human retina using liquid chromatography/tandem mass spectrometry. *J. Chromatogr. A.* **1217(49)**, 7738-7748 (2010).
- Shen, Q., Wang, Y., Gong, L., *et al.* Shotgun lipidomics strategy for fast analysis of phospholipids in fisheries waste and its potential in species differentiation. *J. Agric. Food Chem.* **60(37)**, 9384-9393 (2012).

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, 08/31/2022

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