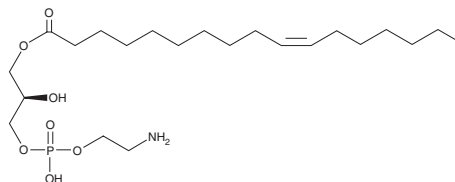


PRODUCT INFORMATION

1-10(Z)-Heptadecenoyl-2-hydroxy-*sn*-glycero-3-PE

Item No. 35672

CAS Registry No.: 1246298-09-8
Formal Name: 10z-heptadecenoic acid, (2R)-3-[[[2-aminoethoxy]hydroxyphosphinyl]oxy]-2-hydroxypropyl ester
Synonyms: LPE(17:1), 17:1 Lyso PE, PE(17:1/0:0)
MF: C₂₂H₄₄NO₇P
FW: 465.6
Purity: ≥98% (9:1 mixture of 1-acyl and 2-acyl)
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

1-10(Z)-Heptadecenoyl-2-hydroxy-*sn*-glycero-3-PE is supplied as a solid. A stock solution may be made by dissolving the 1-10(Z)-heptadecenoyl-2-hydroxy-*sn*-glycero-3-PE in the solvent of choice, which should be purged with an inert gas. 1-10(Z)-Heptadecenoyl-2-hydroxy-*sn*-glycero-3-PE is slightly soluble in chloroform.

Description

1-10(Z)-Heptadecenoyl-2-hydroxy-*sn*-glycero-3-PE is a lysophospholipid. Extracellular levels of 1-10(Z)-heptadecenoyl-2-hydroxy-*sn*-glycero-3-PE in *M. tuberculosis* culture supernatant increase in response to *S. miltiorrhiza* antibacterial tanshinones.¹ It has been used as an internal standard for the quantification of lysophosphoethanolamines.^{2,3}

References

1. Sieniawska, E., Sawicki, R., Marchev, A.S., et al. Tanshinones from *Salvia miltiorrhiza* inhibit *Mycobacterium tuberculosis* via disruption of the cell envelope surface and oxidative stress. *Food Chem. Toxicol.* **156**, 112405 (2021).
2. Medved, J., Knott, B.M., Tarrah, S.N., et al. The lysophospholipid-binding molecule CD1D is not required for the alloimmunization response to fresh or stored RBCs in mice despite RBC storage driving alterations in lysophospholipids. *Transfusion* **61**(7), 2169-2178 (2021).
3. Li, M., Feng, B., Liang, Y., et al. Lipid profiling of human plasma from peritoneal dialysis patients using an improved 2D (NP/RP) LC-QToF MS method. *Anal. Bioanal. Chem.* **405**(21), 6629-6638 (2013).

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, 03/11/2026

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