

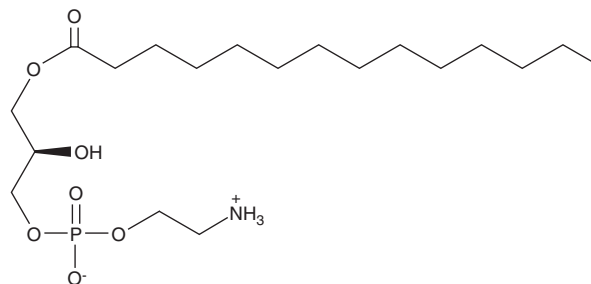
# PRODUCT INFORMATION



## 1-Myristoyl-2-hydroxy-*sn*-glycero-3-PE

Item No. 25591

**CAS Registry No.:** 123060-40-2  
**Formal Name:** tetradecanoic acid, (2R)-3-[[[(2-aminoethoxy)hydroxyphosphinyl]oxy]-2-hydroxypropyl ester  
**Synonyms:** 14:0 Lyso-PE, 1-Myristoyl-2-hydroxy-*sn*-glycero-3-Phosphoethanolamine  
**MF:** C<sub>19</sub>H<sub>40</sub>NO<sub>7</sub>P  
**FW:** 425.5  
**Purity:** ≥98%  
**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-Myristoyl-2-hydroxy-*sn*-glycero-3-PE is supplied as a solid. A stock solution may be made by dissolving the 1-myristoyl-2-hydroxy-*sn*-glycero-3-PE in the solvent of choice. 1-Myristoyl-2-hydroxy-*sn*-glycero-3-PE is soluble in organic solvents such as DMSO, dimethyl formamide, and chloroform, which should be purged with an inert gas.

### Description

1-Myristoyl-2-hydroxy-*sn*-glycero-3-PE is a naturally occurring lysophospholipid.<sup>1</sup> It induces transient increases in intracellular calcium in PC12 cells.<sup>2</sup> Serum levels of 1-myristoyl-2-hydroxy-*sn*-glycero-3-PE are elevated in patients with malignant breast cancer compared to healthy controls.<sup>1</sup>

### References

1. More, T.H., Bagadi, M., RoyChoudhury, S., *et al.* Comprehensive quantitative lipidomic approach to investigate serum phospholipid alterations in breast cancer. *Metabolomics* **13:3** (2017).
2. Lee, J.M., Park, S.J., and Im, D.S. Lysophosphatidylethanolamine increases intracellular Ca<sup>2+</sup> through LPA<sub>1</sub> in PC-12 neuronal cells. *Biochem. Biophys. Res. Commun.* **461(2)**, 378-382 (2015).

#### 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, 12/08/2022

#### 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