

# PRODUCT INFORMATION



## 1,2-Dipalmitoyl-*sn*-glycero-3-N-methyl-PE

Item No. 26010

CAS Registry No.: 3930-13-0

Formal Name: hexadecanoic acid, 1,1'-[(1R)-1-(3-hydroxy-3-oxido-2,4-dioxo-7-aza-3-phosphaoct-1-yl)-1,2-ethanediyl] ester

Synonyms: 1,2-Dipalmitoyl-*sn*-glycero-3-N-methyl-phosphoethanolamine, 16:0 Monomethyl PE, 1,2-NMeDPPE

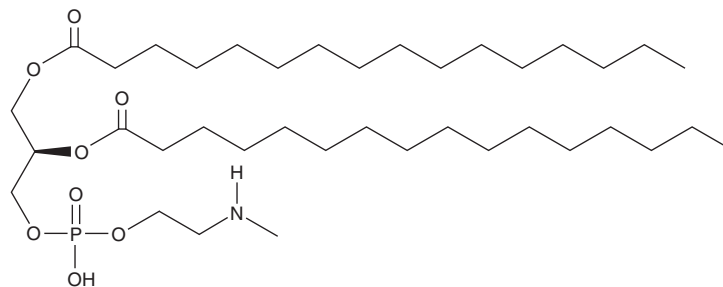
MF: C<sub>38</sub>H<sub>76</sub>NO<sub>8</sub>P

FW: 706.0

Supplied as: A solid

Storage: -20°C

Stability: ≥2 years



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

### Laboratory Procedures

1,2-Dipalmitoyl-*sn*-glycero-3-N-methyl-PE is supplied as a solid. A stock solution may be made by dissolving the 1,2-dipalmitoyl-*sn*-glycero-3-N-methyl-PE in water. We do not recommend storing the aqueous solution for more than one day.

### Description

1,2-Dipalmitoyl-*sn*-glycero-3-N-methyl-PE is a form of 1,2-dipalmitoyl-*sn*-glycero-3-PE (1,2-DPPE; Item No. 15092) that contains a methyl group on the *sn*-3 moiety. The methylated headgroup of 1,2-dipalmitoyl-*sn*-glycero-3-N-methyl-PE does not extend as far into the aqueous subphase as the amine headgroup of 1,2-DPPE when incorporated into a water surface monolayer.<sup>1</sup> It also lowers the phase transition temperature compared with 1,2-DPPE.<sup>2</sup>

### References

1. Dyck, M., Krüger, P., and Lösche, M. Headgroup organization and hydration of methylated phosphatidylethanolamines in Langmuir monolayers. *Phys. Chem. Chem. Phys.* **7**(1), 150-156 (2005).
2. Chowdhry, B.Z. and Dalziel, A.W. Phase transition properties of 1,2- and 1,3-diacylphosphatidylethanolamines with modified head groups. *Biochemistry* **24**(15), 4109-4117 (1985).

#### 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, 08/31/2018

#### 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