

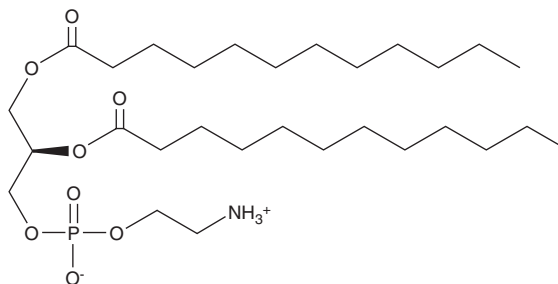
PRODUCT INFORMATION



1,2-Dilauroyl-*sn*-glycero-3-PE

Item No. 15089

CAS Registry No.: 59752-57-7
Formal Name: 1,2-dilauroyl-*sn*-glycero-3-phosphatidylethanolamine
Synonyms: 1,2-Dilauroyl-*sn*-glycero-3-Phosphoethanolamine, DLPE, 1,2-DLPE
MF: C₂₉H₅₈NO₈P
FW: 579.8
Purity: ≥98%
Supplied as: A crystalline 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

1,2-Dilauroyl-*sn*-glycero-3-PE (DLPE) is supplied as a crystalline solid. A stock solution may be made by dissolving the DLPE in the solvent of choice, which should be purged with an inert gas. DLPE is soluble in the organic solvent chloroform at a concentration of approximately 3 mg/ml.

Description

DLPE is a phospholipid containing the medium-chain (12:0) lauric acid inserted at the *sn*-1 and *sn*-2 positions.

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, 12/01/2022

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