

PRODUCT INFORMATION



N-Palmitoyl-DPPE (triethylammonium salt)

Item No. 39195

Formal Name: hexadecanoic acid, (1R)-1-(3-hydroxy-3-oxido-8-oxo-2,4-dioxo-7-aza-3-phosphatricos-1-yl)-1,2-ethanediyl ester, monotriethylammonium salt

Synonyms: N-Palmitoyl-1,2-Dipalmitoyl-*sn*-glycero-3-Phosphoethanolamine, N-Hexadecanoyl-1,2-Dihexadecanoyl-*sn*-glycero-3-Phosphoethanolamine, NAPE(16:0/16:0/16:0)

MF: $C_{53}H_{103}NO_9P \cdot C_6H_{16}N$

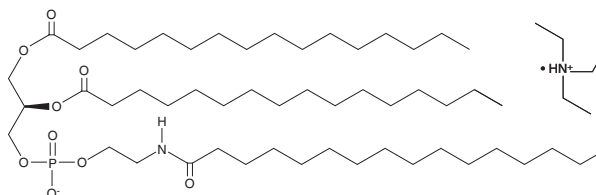
FW: 1,031.6

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

N-Palmitoyl-DPPE (triethylammonium salt) is supplied as a solid. A stock solution may be made by dissolving the N-palmitoyl-DPPE (triethylammonium salt) in the solvent of choice, which should be purged with an inert gas. N-Palmitoyl-DPPE (triethylammonium salt) is soluble in the organic solvent chloroform at a concentration of approximately 5 mg/ml.

Description

N-Palmitoyl-DPPE is an N-acyl-phosphatidylethanolamine (NAPE).¹

Reference

1. Dosoky, N.S., Guo, L., Chen, Z., *et al.* Dietary fatty acids control the species of N-acyl-phosphatidylethanolamines synthesized by therapeutically modified bacteria in the intestinal tract. *ACS Infect. Dis.* **4**(1), 3-13 (2018).

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, 09/17/2024

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