

PRODUCT INFORMATION



1,2-Dipalmitoyl-3-dimethylammonium-propane

Item No. 38311

CAS Registry No.: 96326-74-8

Formal Name: hexadecanoic acid,
1,1'-[1-[(dimethylamino)methyl]-
1,2-ethanediyl] ester

Synonyms: 16:0 DAP, DPDAP

MF: $C_{37}H_{73}NO_4$

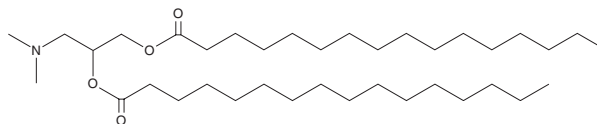
FW: 596.0

Purity: $\geq 95\%$

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

1,2-Dipalmitoyl-3-dimethylammonium-propane (DPDAP) is supplied as a solid. A stock solution may be made by dissolving the DPDAP in the solvent of choice, which should be purged with an inert gas. DPDAP is soluble in the organic solvent chloroform.

Description

DPDAP is an ionizable cationic lipid that has been used in the generation of liposomes for the delivery of miRNA.¹ Liposomes containing DPDAP and encapsulating miR-21, miR-324, and scramble miRNA increase axonogenesis of mouse trigeminal ganglion neurons.

Reference

1. Amit, M., Takahashi, H., Dragomir, M.P., *et al.* Loss of p53 drives neuron reprogramming in head and neck cancer. *Nature* **578(7795)**, 449-454 (2020).

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, 05/05/2023

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