

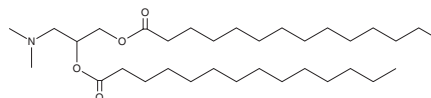
PRODUCT INFORMATION



1,2-Dimyristoyl-3-dimethylammonium-propane

Item No. 38310

CAS Registry No.: 72719-84-7
Formal Name: tetradecanoic acid, 1,1'-[1-[(dimethylamino)methyl]-1,2-ethanediyl] ester
Synonyms: 14:0 DAP, DMDAP
MF: $C_{33}H_{65}NO_4$
FW: 539.9
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-Dimyristoyl-3-dimethylammonium-propane is supplied as a solid. A stock solution may be made by dissolving the 1,2-dimyristoyl-3-dimethylammonium-propane in the solvent of choice, which should be purged with an inert gas. 1,2-Dimyristoyl-3-dimethylammonium-propane is soluble in chloroform.

Description

1,2-Dimyristoyl-3-dimethylammonium-propane is an ionizable cationic lipid that has been used in the generation of liposomes.¹

Reference

1. Nomura, F., Nagata, M., Inaba, T., et al. Capabilities of liposomes for topological transformation. *Chemphyschem* **3**(7), 571-574 (2001).

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