

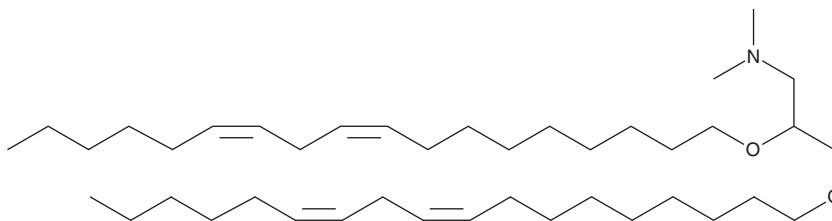
PRODUCT INFORMATION



DLin-DMA

Item No. 36701

CAS Registry No.: 871258-12-7
Formal Name: N,N-dimethyl-2,3-bis[(9Z,12Z)-9,12-octadecadien-1-yloxy]-1-propanamine
Synonym: 1,2-Dilinoleyloxy-N,N-dimethyl-3-aminopropane
MF: C₄₁H₇₇NO₂
FW: 616.1
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

DLin-DMA is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol and dimethyl formamide purged with an inert gas can be used. The solubility of DLin-DMA in these solvents is approximately 10 and 2 mg/ml, respectively.

Description

DLin-DMA is an ionizable cationic amino lipid ($pK_a = 6.7$) that has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs) for the delivery of nucleic acids^{1,2}

References

1. Heyes, J., Palmer, L., Bremner, K., *et al.* Cationic lipid saturation influences intracellular delivery of encapsulated nucleic acids. *J. Control Release* **107(2)**, 276-287 (2005).
2. Kuboyama, T., Yagi, K., Naoi, T., *et al.* Simplifying the chemical structure of cationic lipids for siRNA-lipid nanoparticles. *ACS Med. Chem. Lett.* **10(5)**, 749-753 (2019).

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, 11/08/2022

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