

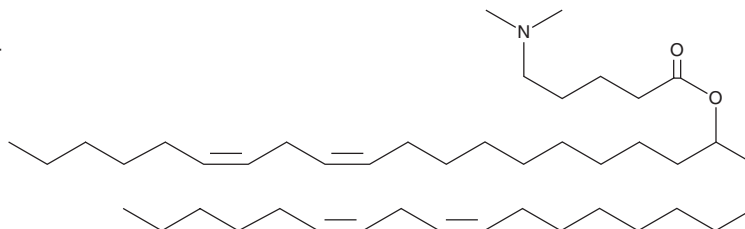
PRODUCT INFORMATION



DLin-MC4-DMA

Item No. 39732

CAS Registry No.: 1226909-66-5
Formal Name: 5-(dimethylamino)-pentanoic acid,
(10Z,13Z)-1-(9Z,12Z)-9,12-octadecadien-
1-yl-10,13-nonadecadien-1-yl ester
Synonym: MC4
MF: C₄₄H₈₁NO₂
FW: 656.1
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-MC4-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. DLin-MC4-DMA is soluble in chloroform and methanol.

Description

DLin-MC4-DMA is an ionizable cationic lipid ($pK_a = 6.93$).¹ Lipid nanoparticles (LNPs) containing DLin-MC4-DMA and encapsulating siRNA for the gene encoding mouse coagulation factor VII (F7) reduce its gene expression levels in mice.

Reference

1. Jayaraman, M., Ansell, S.M., Mui, B.L., *et al.* Maximizing the potency of siRNA lipid nanoparticles for hepatic gene silencing in vivo. *Angew. Chem. Int. Ed. Engl.* **51**(34), 8529-8533 (2012).

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, 11/02/2023

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