

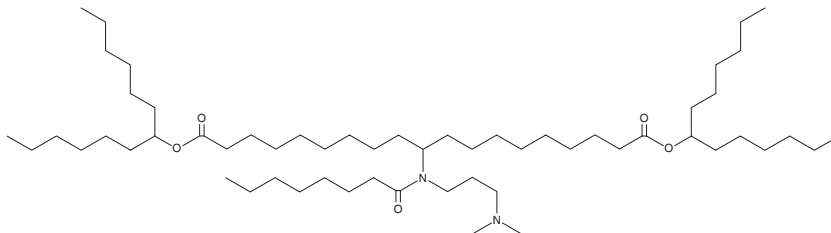
PRODUCT INFORMATION



Lipid 2-10

Item No. 40811

CAS Registry No.: 2055939-68-7
Formal Name: 10-[[3-(dimethylamino)propyl](1-oxooctyl)amino]-nonadecanedioic acid, 1,19-bis(1-hexylheptyl) ester
MF: C₅₈H₁₁₄N₂O₅
FW: 919.5
Purity: ≥98%
Supplied as: A 5 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Lipid 2-10 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. Lipid 2-10 is sparingly soluble (1-10 mg/ml) in chloroform and methanol.

Description

Lipid 2-10 is an ionizable cationic lipid that has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs).¹

Reference

1. Hope, M.J., Mui, B., Lin, P.J.C., *et al.* Lipid nanoparticle formulations. *Acuitas Therapeutics, Inc.* US 11,712,481 B2, (2023).

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, 06/25/2026

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