

PRODUCT INFORMATION



CP-LC-1254

Item No. 41928

CAS Registry No.: 3040859-52-4

Formal Name: 2-[[[2-(dimethylamino)ethyl]amino]carbonyl]-4-[[3-[(2-hexyldecyl)oxy]-3-oxopropyl]thio]-butanoic acid, 2-hexyldecyl ester

MF: C₄₄H₈₆N₂O₅S

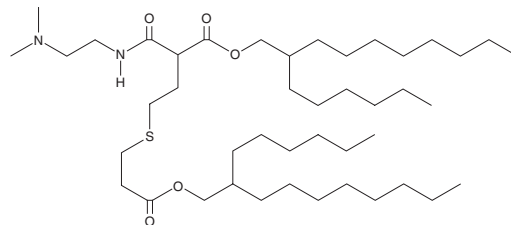
FW: 755.2

Purity: ≥98%

Supplied as: A 10 mg/ml solution in ethanol

Storage: -20°C

Stability: ≥1 year



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

Description

CP-LC-1254 is an ionizable cationic lipid that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of circular RNA (circRNA) *in vivo*.¹ Intramuscular administration of LNPs containing CP-LC-1254 and encapsulating circRNA encoding a luciferase reporter induces luciferase expression in mice.

Reference

1. Broset, E., Larraga, A., Lampaya, V., et al. A complete approach for circRNA therapeutics from purification to lyophilized delivery using novel ionizable lipids. *Int. J. Mol. Sci.* **26(11)**, 5138 (2025).

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, 11/21/2025

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