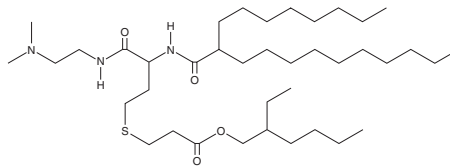


PRODUCT INFORMATION



CP-LC-0867
Item No. 41926

CAS Registry No.: 3040858-96-3
Formal Name: 3-[[4-[[2-(dimethylamino)ethyl]amino]-3-[[2-octyl-1-oxododecyl]amino]-4-oxobutyl]thio]-propanoic acid, 2-ethylhexyl ester
MF: C₃₉H₇₇N₃O₄S
FW: 684.1
Purity: ≥98%
Supplied as: A 50 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-0867 is an ionizable cationic lipid that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of circular RNA (circRNA) and mRNA *in vivo*.^{1,2} Intramuscular administration of LNPs containing CP-LC-0867 and encapsulating circRNA encoding a luciferase reporter induce the production of luciferase protein for at least 14 days in mice.² LNP-based vaccines containing CP-LC-0867 and encapsulating mRNA encoding the severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) spike protein increase B- and T cell responses in mice with an increase in serum IgG titers and splenocyte IFN-γ levels, respectively, after six weeks.¹ CP-LC-0867-containing LNP-based mRNA vaccines increase splenocyte IFN-γ levels in mice to a greater extent than vaccines containing CP-LC-0431 or CP-LC-0474.

References

1. Mata, E., Broset, E., Matute, C., *et al.* Preserved efficacy of lyophilized SARS-CoV-2 mRNA vaccine incorporating novel ionizable lipids after one year at 25 °C. *NPJ Vaccines* **10**(1), 135 (2025).
2. 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, 05/04/2026

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