

PRODUCT INFORMATION



CP-LC-1447
Item No. 41922

CAS Registry No.: 3112948-20-3
Formal Name: 14-[[[2-(1H-imidazol-1-yl)ethyl]amino]carbonyl]-7-[[[3-(octyloxy)-3-oxopropyl]thio]-12,20-dioxo-21-oxa-4,17-dithia-13-azatetratriacontanoic acid, octyl ester

MF: C₅₅H₁₀₀N₄O₈S₃

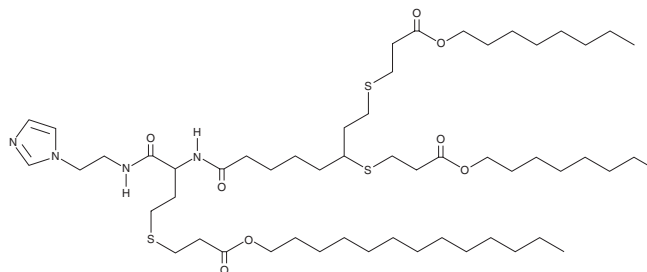
FW: 1,041.6

Purity: ≥95%

Supplied as: A 10 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.

Description

Lipid CP-LC-1447 is an ionizable cationic amino lipid derived from homocysteine, a naturally occurring amino acid. This lipid has demonstrated an efficient delivery and high protein expression of different kinds of RNA (mRNA, cRNA and saRNA) *in vivo*. Moreover, using a typical LNP formulation (IL/Ch/DOPE/PEG: 50/38.5/10/1.5), it has shown to induce specific spleen targeting much more efficiently than other commercial alternatives through IV injection, with no signs of toxicity.

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, 07/02/2026

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