

PRODUCT INFORMATION



CP-LC-0729

Item No. 41929

CAS Registry No.: 3040858-60-1

Formal Name: 3-[[4-[[2-(dimethylamino)ethyl]amino]-3-[(2-hexyl-1-oxodecyl)amino]-4-oxobutyl]thio]-propanoic acid, 2-hexyldecyl ester

Synonym: A4B2C8

MF: C₄₃H₈₅N₃O₄S

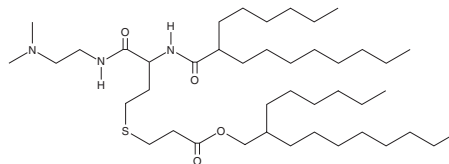
FW: 740.2

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-0729 is an ionizable cationic amino lipid (apparent $pK_a = 6.78$) that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vivo*.¹ Intravenous administration of LNPs containing CP-LC-0729 and encapsulating mRNA encoding luciferase selectively accumulate in the lungs and, to a lesser extent, in the spleen over the kidneys, liver, heart, or intestines in mice.

Reference

1. Peña, Á., Heredero, J., Blandín, B., *et al.* Multicomponent thiolactone-based ionizable lipid screening platform for efficient and tunable mRNA delivery to the lungs. *Commun. Chem.* **8**(1), 116 (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, 04/21/2026

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