

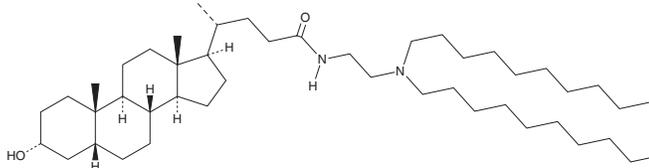
PRODUCT INFORMATION



LC10

Item No. 37000

CAS Registry No.: 2768690-22-6
Formal Name: (3 α ,5 β)-N-[2-(didecylamino)ethyl]-3-hydroxy-cholan-24-amide
MF: C₄₆H₈₆N₂O₂ • HCl
FW: 735.7
Purity: ≥98%
Supplied as: A 100 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

LC10 is a lipidoid derivative of the secondary bile acid lithocholic acid (Item No. 20253) that contains two 10-carbon alkyl chains.¹ It has been used in the generation of self-assembling skin-permeable lipid nanoparticles (LNPs). LC10 LNPs inhibit the proliferation of normal human epidermal keratinocytes (NHEKs) when used at a concentration of 100 nM and reduce the levels of TNF- α , IL-17, and IL-23 in human peripheral blood mononuclear cells (PBMCs) at 50, 100, and 200 nM. Topical application of LC10 LNPs (10 mg/kg) reduces psoriatic-like plaques and reverses weight loss induced by imiquimod (Item No. 14956) in a mouse model of psoriasis.

Reference

1. Rachamalla, H.K., Voshavar, C., Arjunan, P., *et al.* Skin-permeable nano-lithocholic lipidoid efficiently alleviates psoriasis-like chronic skin inflammations. *ACS Appl. Mater. Interfaces* **14(13)**, 14859-14870 (2022).

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, 12/01/2025

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