

# PRODUCT INFORMATION



## H1L1A1B3

Item No. 41199

**Formal Name:** 7-(cyclohexylamino)-6-(N-(3-(dimethylamino)propyl)-4-methylnonanamido)-7-oxoheptyl dodecanoate

**MF:** C<sub>40</sub>H<sub>77</sub>N<sub>3</sub>O

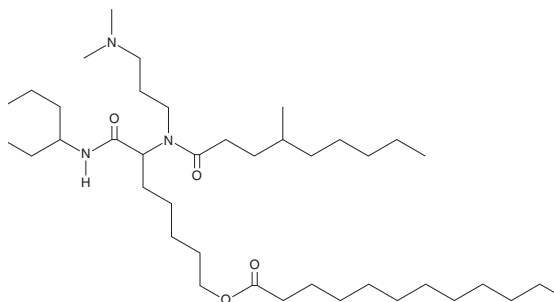
**FW:** 664.1

**Purity:** ≥95%

**Supplied as:** A 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

H1L1A1B3 is an ionizable cationic lipid that has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs) for the delivery of circular RNA (circRNA) *in vitro* and *in vivo*.<sup>1</sup> LNPs containing H1L1A1B3 and encapsulating circRNA encoding IL-12 induce IL-12 production four times more effectively than LNPs containing ALC-0315 (Item No. 34337) and encapsulating the same circRNA in LLC1 murine Lewis lung carcinoma cells. Administration of LNPs containing H1L1A1B3 and encapsulating circRNA encoding IL-12 reduces tumor growth in an LLC1 murine Lewis lung carcinoma model in combination with an antibody against programmed cell death ligand 1 (PD-L1).

### Reference

1. Xu, S., Xu, Y., Solek, N.C., *et al.* Tumor-tailored ionizable lipid nanoparticles facilitate IL-12 circular RNA delivery for enhanced lung cancer immunotherapy. *Adv. Mater.* **36(29)**, e2400307 (2024).

#### 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, 09/10/2024

#### 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