

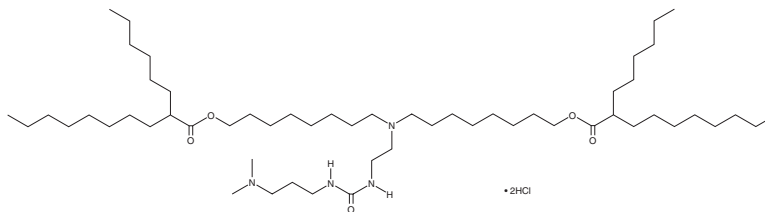
PRODUCT INFORMATION



Lipid 35 (hydrochloride)

Item No. 43127

Formal Name: 11-(8-((2-hexyldecanoyl)oxy)octyl)-2-methyl-7-oxo-2,6,8,11-tetraazanonadecan-19-yl 2-hexyldecanoate, dihydrochloride
MF: C₅₆H₁₁₂N₄O₅ • 2HCl
FW: 994.5
Purity: ≥95%
Supplied as: A 10 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Lipid 35 is an ionizable cationic lipid ($pK_a = 7.6$) that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vitro* and *in vivo*.¹ LNPs containing lipid 35 selectively accumulate in the lungs, including in pulmonary epithelial, endothelial, myeloid, and dendritic cells, and spleen in mice. LNPs containing lipid 35 and encapsulating mRNA encoding *Pseudomonas* exotoxin A decrease the number of lung metastases and increase survival in a B16/F10.9 murine melanoma model.

Reference

1. Naidu, G.S., Rampado, R., Sharma, P., *et al.* Ionizable lipids with optimized linkers enable lung-specific, lipid nanoparticle-mediated mRNA delivery for treatment of metastatic lung tumors. *ACS Nano* **19**(6), 6571-6587 (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, 06/24/2026

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