

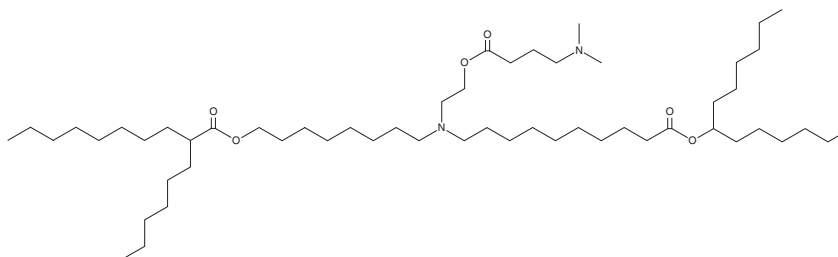
PRODUCT INFORMATION



Lipid 23

Item No. 38862

CAS Registry No.: 3117189-61-1
Formal Name: 8-((2-((4-(dimethylamino)butanoyl)oxy)ethyl)(10-oxo-10-(tridecan-7-yloxy)decyl)amino)octyl 2-hexyldecanoate
MF: C₅₅H₁₀₈N₂O₆
FW: 893.5
Purity: ≥95%
Supplied as: A 25 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 23 is an ionizable cationic amino lipid ($pK_a = 5.7$) that has been used with other lipids in the formulation of lipid nanoparticles (LNPs).¹ LNPs containing lipid 23 and encapsulating an mRNA reporter accumulate specifically in mouse liver after intravenous administration.

Reference

1. Naidu, G.S., Yong, S.-B., Ramishetti, S., *et al.* A combinatorial library of lipid nanoparticles for cell type-specific mRNA delivery. *Adv. Sci. (Weinh)* **10(19)**, e2301929 (2023).

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, 05/04/2026

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