

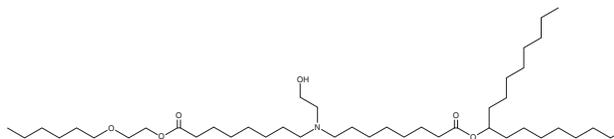
PRODUCT INFORMATION



Lipid ND-O1

Item No. 43010

CAS Registry No.: 3096665-79-8
Formal Name: heptadecan-9-yl 8-((8-(2-(hexyloxy)ethoxy)-8-oxooctyl)(2-hydroxyethyl)amino)octanoate
Synonym: SM-86 Analog-2
MF: C₄₃H₈₅NO₆
FW: 712.2
Purity: ≥95%
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Description

Lipid ND-O1 is an ether-containing ionizable cationic lipid ($pK_a = 6.6$) that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of siRNA *in vitro* and *in vivo*.¹ LNPs containing ND-O1 and encapsulating Cy5-labeled siRNA selectively accumulate in the liver over the spleen, heart, kidney, and lung in mice. LNPs containing ND-O1 and encapsulating siRNA targeting heat shock protein 47 (Hsp47) decrease hepatic collagen, TNF- α , IL-1 β , TGF- β , and IL-6 levels in a mouse model of hepatic fibrosis induced by carbon tetrachloride (CCl₄).

Reference

1. Chu, R., Kong, J., Gao, Q., *et al.* Ether bond-modified lipid nanoparticles for enhancing the treatment effect of hepatic fibrosis. *Int. J. Pharm.* **671**, 125192 (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, 10/13/2025

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