

PRODUCT INFORMATION



306-O12B-3

Item No. 37096

CAS Registry No.: 2982842-90-8
Formal Name: bis(2-(octyldisulfaneyl)ethyl)
3,3'-((4-methyl-11-oxo-12-oxa-
15,16-dithia-4,8-diazatetracosyl)
azanediyldipropionate

MF: C₄₆H₉₁N₃O₆S₆

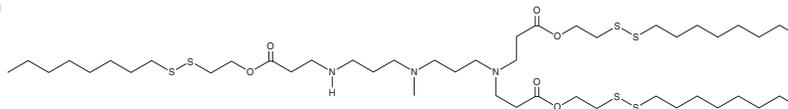
FW: 974.6

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

306-O12B-3 is an ionizable cationic lipidoid.¹ It has been used in the generation of lipid nanoparticles (LNPs) for the delivery of antisense oligonucleotides (ASOs). LNPs containing 306-O12B-3 localize to the liver after intravenous administration in mice. LNPs containing 306-O12B-3 in combination with other lipids encapsulating ASOs targeting the gene encoding proprotein convertase subtilisin kexin type 9 (PCSK9) decrease hepatic PCSK9 expression in mice. Intravenous administration of LNPs containing 306-O12B-3 and the cationic lipidoid NT1-O14B (Item No. 37095) encapsulating ASOs against tau decreases tau expression and levels in mouse brain.²

References

1. Yang, L., Ma, F., Liu, F., *et al.* Efficient delivery of antisense oligonucleotides using bioreducible lipid nanoparticles *in vitro* and *in vivo*. *Mol. Ther. Nucleic Acids* **19**, 1357-1367 (2020).
2. Ma, F., Yang, L., Sun, Z., *et al.* Neurotransmitter-derived lipidoids (NT-lipidoids) for enhanced brain delivery through intravenous injection. *Sci. Adv.* **6**(30), eabb4429 (2020).

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, 11/06/2025

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