

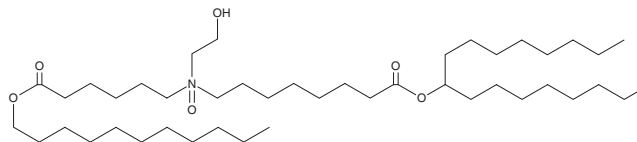
PRODUCT INFORMATION



SM-102 N-oxide

Item No. 39636

CAS Registry No.: 2824195-50-6
Formal Name: 8-(heptadecan-9-yloxy)-N-(2-hydroxyethyl)-8-oxo-N-(6-oxo-6-(undecyloxy)hexyl)octan-1-amine oxide
Synonyms: CAY10779 N-oxide, Lipid H N-oxide, LNP-102 N-oxide
MF: C₄₄H₈₇NO₆
FW: 726.2
Purity: ≥98%
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

SM-102 N-oxide is potential impurity in commercial preparations of SM-102.¹

References

1. Packer, M., Gyawali, D., Hua, S., et al. Lipid nanoparticle compositions and methods of formulating the same. *EP 4 046 629 A1* (2022).

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, 10/10/2023

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