

PRODUCT INFORMATION



SM-102

Item No. 33474

CAS Registry No.: 2089251-47-6
Formal Name: 8-[(2-hydroxyethyl)[6-oxo-6-(undecyloxy)hexyl]amino]-octanoic acid, 1-octylnonyl ester

Synonyms: CAY10779, Lipid H, LNP-102

MF: C₄₄H₈₇NO₅

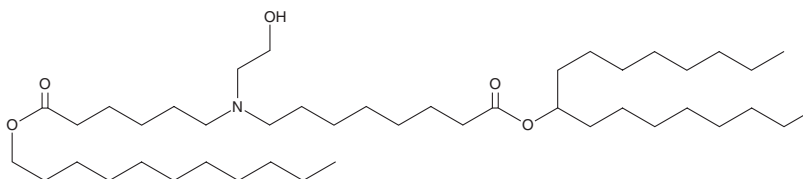
FW: 710.2

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 is an ionizable cationic amino lipid that has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs).¹ Administration of luciferase mRNA in SM-102-containing LNPs induces hepatic luciferase expression in mice. Formulations containing SM-102 have been used in the development of LNPs for delivery of mRNA-based vaccines.

Reference

1. Sabnis, S., Kumarasinghe, E.S., Salerno, T., *et al.* A novel amino lipid series for mRNA delivery: Improved endosomal escape and sustained pharmacology and safety in non-human primates. *Mol. Ther.* **26(6)**, 1509-1519 (2018).

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, 03/29/2023

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