

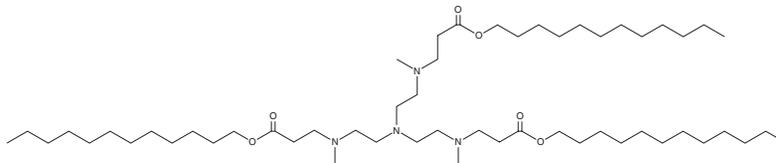
PRODUCT INFORMATION



SIL Lipid

Item No. 40801

CAS Registry No.: 1566559-79-2
Formal Name: N,N',N''-(nitrilotri-2,1-ethanediyl) tris[N-methyl-β-alanine, 1,1',1''-tridodecyl ester]
MF: C₅₄H₁₀₈N₄O₆
FW: 909.5
Purity: ≥98%
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

SIL lipid is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice, which should be purged with an inert gas. SIL lipid is sparingly soluble (1-10 mg/ml) in DMSO.

Description

SIL lipid is an ionizable cationic lipid that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of siRNA *in vitro*.¹ LNPs containing SIL lipid and encapsulating siRNA targeting mRNA encoding the purinergic P2X₇ receptor decrease protein levels of the purinergic P2X₇ receptor, reduce migration in a wound healing assay, and induce apoptosis and necrosis in 4T1 breast cancer cells.

Reference

1. Kiaie, S.H., Zangi, A.R., Sheibani, M., *et al.* Novel synthesized ionizable lipid for LNP-mediated P2X₇siRNA to inhibit migration and induce apoptosis of breast cancer cells. *Purinergic Signal* (2024).

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, 11/18/2025

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