

PRODUCT INFORMATION



244cis

Item No. 39284

CAS Registry No.: 2956402-64-3
Formal Name: 4-[2-[bis[9-[(2Z)-2-nonen-1-yloxy]-9-oxononyl]amino]ethyl]-1-piperazinenonanoic acid, (2Z)-2-nonen-1-yl ester

MF: C₆₀H₁₁₁N₃O₆

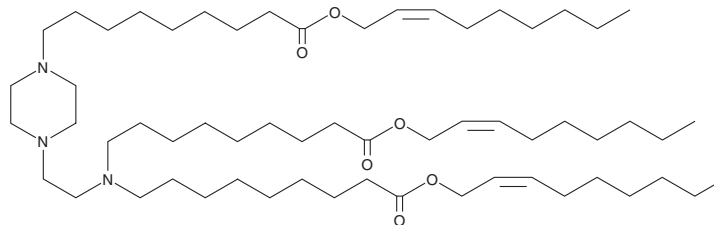
FW: 970.6

Purity: ≥95%

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

244cis is an ionizable cationic lipid that has been used in the generation of lipid nanoparticles (LNPs).¹ LNPs containing 244cis and encapsulating an mRNA reporter specifically accumulate in the lungs and induce lower serum chemokine (C-C motif) ligand 2 (CCL2) levels than LNPs containing SM-102 (Item No. 33474) in mice. Intravenous administration of LNPs containing 244cis and encapsulating mRNA encoding Cre recombinase increases tdTomato fluorescence in lung endothelial, epithelial, and immune cells, as well as fibroblasts, in mice expressing a floxed STOP codon to prevent tdTomato expression.

Reference

1. Kim, M., Jeong, M., Lee, G., *et al.* Novel piperazine-based ionizable lipid nanoparticles allow the repeated dose of mRNA to fibrotic lungs with improved potency and safety. *Bioeng. Transl. Med.* e10556 (2023).

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, 12/11/2023

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