

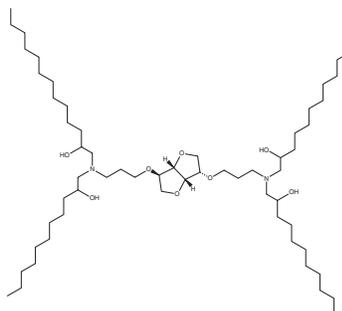
PRODUCT INFORMATION



Lipid DIM1

Item No. 40512

CAS Registry No.: 2820155-36-8
Formal Name: 1,1',1'',1'''-((((3R,3aS,6S,6aS)-hexahydrofuro[3,2-b]furan-3,6-diyl)bis(oxy))bis(propane-3,1-diyl))bis(azanetriyl)tetrakis(dodecan-2-ol)
MF: C₆₀H₁₂₀N₂O₈
FW: 997.6
Purity: ≥95%
Supplied as: A 50 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

Lipid DIM1 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. Lipid DIM1 is sparingly soluble (1-10 mg/ml) in chloroform and slightly soluble (0.1-1 mg/ml) in methanol purged with an inert gas.

Description

Lipid DIM1 is an ionizable cationic lipid that has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs) for mRNA delivery *in vitro*.¹ LNPs containing lipid DIM1 and encapsulating mRNA encoding a luciferase reporter induce luminescence in primary mouse adipose stem cells (ASCs).

Reference

1. Xue, Y., Zhang, Y., Zhong, Y., *et al.* LNP-RNA-engineered adipose stem cells for accelerated diabetic wound healing. *Nat. Commun.* **15**(1), 739 (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