

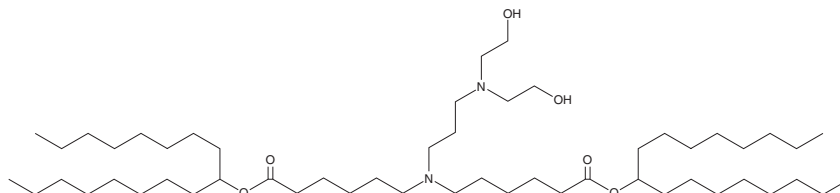
PRODUCT INFORMATION



Lipid U 105

Item No. 41944

CAS Registry No.: 3049129-14-5
Formal Name: di(heptadecan-9-yl) 6,6'-((3-(bis(2-hydroxyethyl)amino)propyl) azanediyl)dihexanoate
MF: C₅₃H₁₀₆N₂O₆
FW: 867.4
Purity: ≥95%
Supplied as: A 10 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Description

Lipid U 105 is an ionizable cationic lipid ($pK_a = 6.65$) that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of circular RNA *in vitro* and *in vivo*.¹ LNPs containing lipid U 105 and encapsulating circular RNA encoding VEGF-A increase the levels of VEGF-A in, and increases the proliferation and migration of, human umbilical vein endothelial cells (HUVECs). Topical administration of the same LNPs increases wound healing in a mouse model of diabetes induced by a high-fat and -sugar diet.

Reference

1. Liu, J., Zhang, Y., Liu, C., *et al.* A single dose of VEGF-A circular RNA sustains in situ long-term expression of protein to accelerate diabetic wound healing. *J. Control Release* **373**, 319-335 (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, 11/21/2025

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