

PRODUCT INFORMATION



C14-306

Item No. 45262

CAS Registry No.: 2153563-36-9

Formal Name: 1,1',1'',1'''-[(methylimino)bis(3,1-propanediyl)nitrido]]tetrakis-2-tetradecanol

MF: $C_{63}H_{131}N_3O_4$

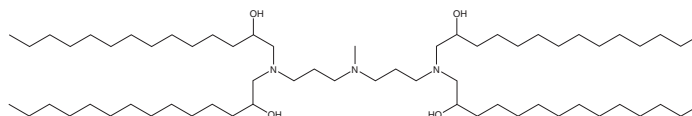
FW: 994.8

Purity: ≥95% (mixture of isomers)

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.

Description

C14-306 is an ionizable cationic lipid that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vitro* and *in vivo*.¹ LNPs containing C14-306 and encapsulating a luciferase mRNA reporter induce luminescence to a greater extent in SH-SY5Y cells than in hCMEC/D3 cells. Intravenous administration of LNPs containing C14-306 accumulate in the mouse brain, liver, and spleen.

Reference

1. Han, E.L., Kim, D., Murray, A.M., *et al.* High-throughput *in vivo* screening identifies structural factors driving mRNA lipid nanoparticle delivery to the brain. *ACS Nano* **20**(4), 3807-3820 (2026).

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, 03/13/2026

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