

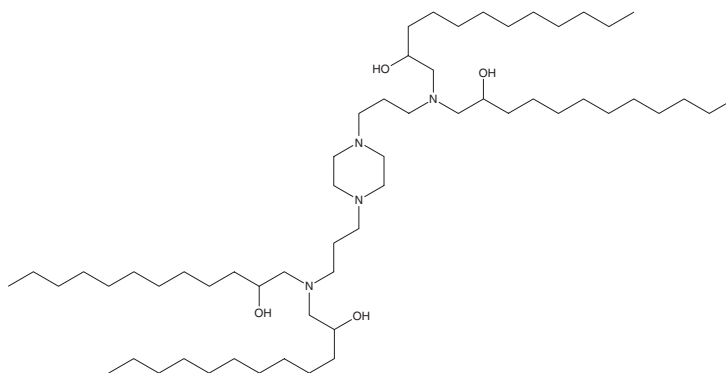
# PRODUCT INFORMATION



**246C10**

Item No. 37907

**CAS Registry No.:** 2635329-26-7  
**Formal Name:** 1,1',1'',1'''-((piperazine-1,4-diylbis(propane-3,1-diyl))bis(azanetriyl))tetrakis(dodecan-2-ol)  
**MF:**  $C_{58}H_{120}N_4O_4$   
**FW:** 937.6  
**Purity:**  $\geq 95\%$   
**Supplied as:** A solution in ethanol  
**Storage:**  $-20^{\circ}\text{C}$   
**Stability:**  $\geq 2$  years



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

## Laboratory Procedures

246C10 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. 246C10 is soluble in the organic solvent DMSO.

## Description

246C10 is an ionizable cationic lipid ( $pK_a = 6.75$ ).<sup>1</sup> It has been used in combination with other lipids in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vitro* and *in vivo*. LNPs containing 246C10 and encapsulating an mRNA reporter selectively accumulate in mouse liver after intravenous administration.

## Reference

1. Kim, M., Jeong, M., Hur, S., *et al.* Engineered ionizable lipid nanoparticles for targeted delivery of RNA therapeutics into different types of cells in the liver. *Sci. Adv.* **7**(9), eabf4398 (2021).

### 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, 03/18/2025

## 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](http://WWW.CAYMANCHEM.COM)