

PRODUCT INFORMATION

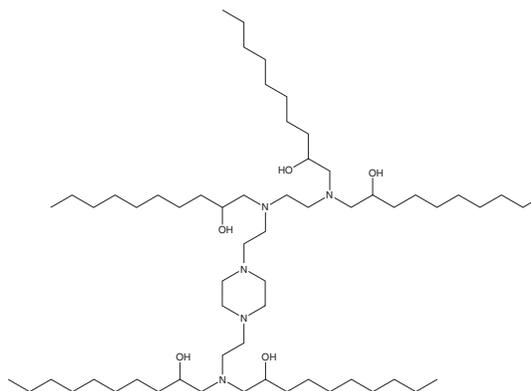


C10-200

Item No. 40863

CAS Registry No.: 3082233-50-6
Formal Name: 1,1'-((2-(4-(2-((2-(bis(2-hydroxydecyl)amino)ethyl)(2-hydroxydecyl)amino)ethyl)piperazin-1-yl)ethyl)azanediyl)bis(decan-2-ol)

MF: C₆₀H₁₂₅N₅O₅
FW: 996.7
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.

Laboratory Procedures

C10-200 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. C10-200 is sparingly soluble (1-10 mg/ml) in chloroform.

Description

C10-200 is an ionizable cationic lipid that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vivo*.¹ LNPs containing C10-200 and encapsulating mRNA encoding a luciferase reporter selectively induce luminescence in the livers of mice to a greater extent than LNPs containing the branched-chain ionizable cationic lipidoid C12-200 (Item No. 36699). LNPs containing C10-200 and encapsulating mRNA encoding erythropoietin (EPO) increase serum levels of EPO in mice.

Reference

1. Mrksich, K., Padilla, M.S., Joseph, R.A., *et al.* Influence of ionizable lipid tail length on lipid nanoparticle delivery of mRNA of varying length. *J. Biomed. Mater. Res. A* (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