

PRODUCT INFORMATION



C9-200

Item No. 40862

Formal Name: 1,1'-((2-(4-(2-((2-(bis(2-hydroxynonyl)amino)ethyl)(2-hydroxynonyl)amino)ethyl)piperazin-1-yl)ethyl)azanediyl)bis(nonan-2-ol)

MF: $C_{55}H_{115}N_5O_5$

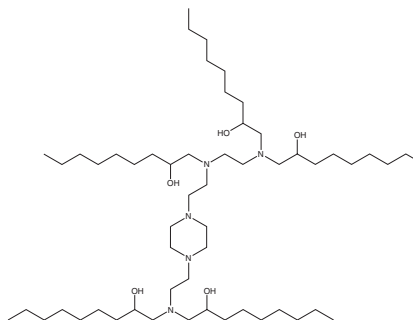
FW: 926.5

Purity: ≥95% (mixture of isomers)

Supplied as: A 10 mg/ml solution in ethanol

Storage: -20°C

Stability: ≥3 years



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

Description

C9-200 is an ionizable cationic lipid that has been used in the formation of lipid nanoparticles (LNPs).¹ LNPs containing C9-200 and encapsulating mRNA encoding erythropoietin (EPO) increase serum EPO levels in mice. LNPs containing C9-200 and encapsulating mRNA encoding the Cas9 nuclease and small-guide RNA (sgRNA) targeting transthyretin (TTR) induce 3-fold higher hepatic insertion and deletion (indel) formation than LNPs containing C12-200 (Item No. 36699) and encapsulating mRNA encoding Cas9 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* **112(9)**, 1494-1505 (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