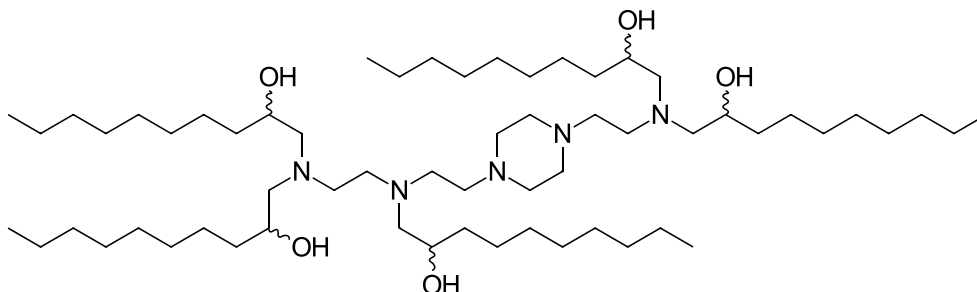


C10-200

Catalog Number: N-1221



Molecular Formula: C₆₀H₁₂₅N₅O₅

MW: 996.66

CAS#: 3082233-50-6

Alternative Names: 1-[N-(2-hydroxydecyl)(2-{N-[2-(4-{2-[bis(2-hydroxydecyl)amino]ethyl}piperazin-1-yl)ethyl]}(2-hydroxydecyl)amino}ethyl)amino]decan-2-ol

Solubility: 50 mg/mL ethanol or methanol, 10 mg/mL chloroform

Storage and Handling: Store dry at -20°C, Stock solution should be stored frozen (-20°C or below).

Background: C10-200 is an ionizable lipid used to generate lipid nanoparticles (LNPs) for delivery of RNA *in vitro* and *in vivo*. C10-200 LNPs formulated with firefly luciferase mRNA showed a 10-fold increase in liver transfection compared with the longer tail analog, C12-200 (Cat# N-1220).

References: Mrksich, K., et al. (2024) "Influence of ionizable lipid tail length on lipid nanoparticle delivery of mRNA of varying length." J. Biomed. Mater. Res. A 112(9), 1494-1505. DOI: 10.1002/jbm.a.37705

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the SDS on request. This product is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.