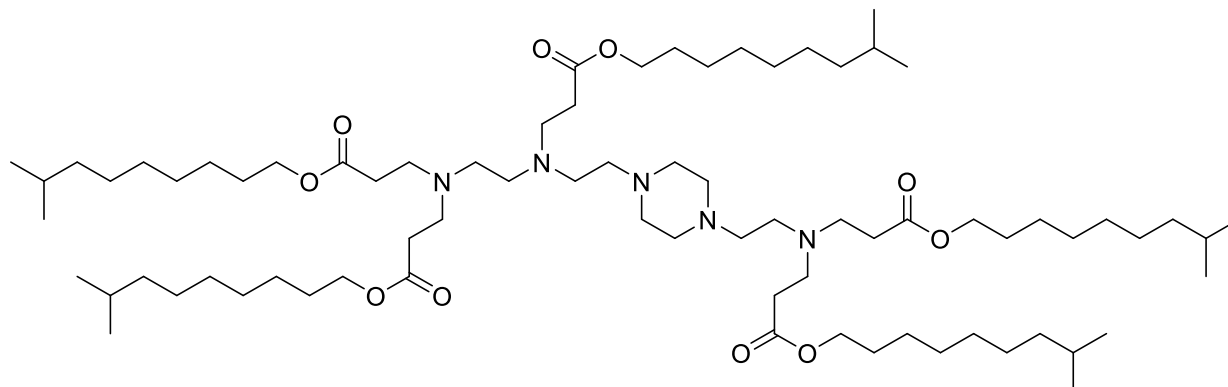


2000i10

Catalog Number: N-1021



Molecular Formula: $C_{75}H_{145}N_5O_{10}$

MW: 1277.01

CAS: 2764986-99-2

Alternative Names: bis(8-methylnonyl) 3,3'-((2-(4-(2-((2-(bis(3-((8-methylnonyl)oxy)-3-oxopropyl)amino)ethyl)(3-((8-methylnonyl)oxy)-3-oxopropyl)amino)ethyl)piperazin-1-yl)ethyl)azanediyl)dipropionate

Solubility: Methanol, ethanol, DMSO.

Storage and Handling: 2000i10 is stable for at least 1 year when stored as oil, protected from moisture, and light at -20°C.

Background: Lipid nanoparticles (LNPs) facilitate the delivery of mRNA. 2000i10 lipid nanoparticles induce protein expression in the liver and the spleen.

References: 1) Jilian R. Melamed et al. "Lipid nanoparticle chemistry determines how nucleoside base modifications alter mRNA delivery." *Journal of Controlled Release* ,2022,341:206-214. doi: <https://doi.org/10.1016/j.jconrel.2021.11.022>

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS 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.

Technical Data Sheet, Rev 1, 09032025 – **For research use only.** Not intended for diagnostic or therapeutic use.