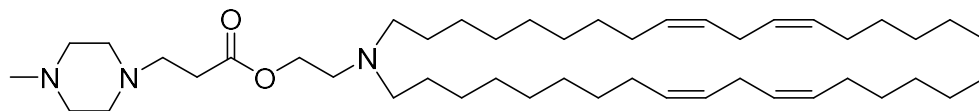


Lipid 10

Catalog number: N-1110



Molecular Formula: C₄₆H₈₅N₃O₂

MW: 712.21

CAS: 2430034-02-7

Alternate Name: 2-(di((9Z,12Z)-octadeca-9,12-dien-1-yl)amino)ethyl 3-(4-methylpiperazin-1-yl)propanoate

Solubility: Methanol, ethanol, DMSO.

Storage and Handling: Lipid 10 is stable for at least 2 years when stored as an oil, protected from moisture, and light at -20°C. Solutions of lipid 10 should be stored at -20°C between uses.

Background: Lipid 10 is an ionizable lipid which efficiently delivers siRNA into leukocytes.

References: 1) Srinivas Ramishetti, et al "A Combinatorial Library of Lipid Nanoparticles for RNA Delivery to Leukocytes" Adv. Mater. 2020,32(12):e1906128. doi: 10.1002/adma.201906128.

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, 5/21/2024 – **For research use only.** Not intended for diagnostic or therapeutic use.