

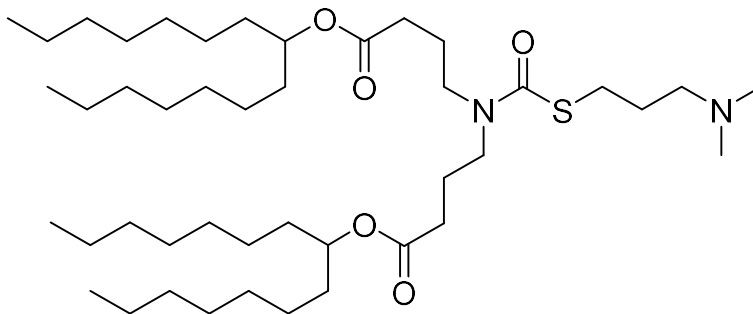
ATX100

Catalog number: N-1100

Molecular Formula: C₄₄H₈₆N₂O₅S

MW: 755.24

CAS#: 2230647-37-5



Alternate Name: Bis(2-(decyldisulfanyl)ethyl) 3,3'-((2-(1H-indol-3-yl)ethyl)azanediyl)dipropionate

Solubility: Methanol, ethanol, DMSO.

Storage and Handling: ATX100 is stable for at least 2 years when stored as an oil, protected from moisture, and light at -20°C. Solutions of ATX100 should be stored at -20°C between uses.

Background: ATX100 is a biodegradable ionizable lipid with improved *in vivo* activities for siRNA delivery via lipid nanoparticles (LNP). The physicochemical parameters of ATX100: pKa 6.316, lipophilicity cLogD 14.03, nanoparticle size 82 nm. The *in vivo* factor VII knockdown efficiency of ATX100 is 97% (0.03mg/kg in mice, i.v.).

References: 1) Kumar Rajappan et al. "Property-Driven Design and Development of Lipids for Efficient Delivery of siRNA" J. Med. Chem. 2020, 63(21):12992-13012. DOI: 10.1021/acs.jmedchem.0c01407

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.