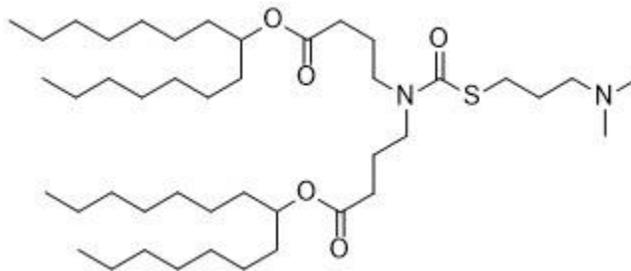


ATX-0126 (ATX100)

Catalog Number: N-1100



Molecular Formula: C₄₄H₈₆N₂O₅S

MW: 755.24

CAS#: 2230647-37-5

Alternative Names: ATX-126, ATX-100, ATX100, Lipid 10p, Bis(2-(decyldisulfanyl)ethyl) 3,3'-((2-(1H-indol-3-yl)ethyl)azanediyl)dipropionate

Solubility: Methanol, ethanol, DMSO.

Storage and Handling: ATX-0126 is stable for at least 2 years when stored as an oil at -20°C, protected from moisture and light. Solutions of ATX-0126 should be maintained at -20°C between uses to ensure stability.

Background: ATX-0126 is a newly developed ionizable, cationic, biodegradable lipid designed for lipid nanoparticle formulations to enhance siRNA delivery. ATX-0126 exhibits: pKa 6.316, cLogD 14.03, average nanoparticle size 82 nm. ATX-0126 induced 97% *in vivo* factor VII knockdown at 0.03 mg/kg (*i.v.*, mice).

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 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.