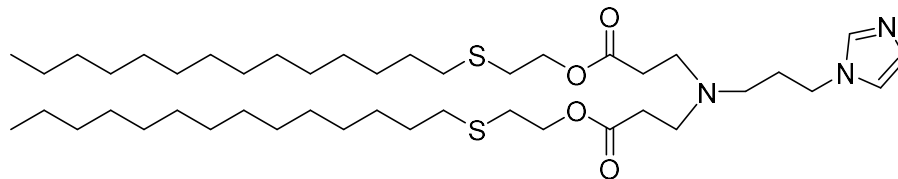


# Echelon Biosciences Inc.

---

## 93-017S

Catalog number: N-1175



**Molecular Formula:** C<sub>44</sub>H<sub>83</sub>N<sub>3</sub>O<sub>4</sub>S<sub>2</sub>

**MW:** 782.29

**CAS:** 2227008-67-3

**Alternative Names:** Bis(2-(tetradecylthio)ethyl 3,3'-((3-(1H-imidazol-1-yl)propyl)azanediyl)dipropionate

**Solubility:** Ethanol, chloroform, Ethyl Acetate

**Storage and Handling:** Store dry at -20 °C or below. Stock solutions should be stored at -20 °C.

**Background:** 93-017S is a lipidoid used for generating lipid nanoparticle to deliver RNA, peptides, and nucleotides. 93-017S LNPs were able to deliver Cre recombinase mRNA in mice. In addition, LNPs were able to deliver the Cre recombinase protein and Cas9: single-guide RNA into cells for genome editing. 93-017S LNPs were used to deliver the STING pathway agonist 2'3'-Cyclic GMP-AMP (cGAMP) when injected into B16F10 tumors in mice which has potential as an *in situ* cancer vaccine.

**References:** 1) Zhao X, Chen J, Qiu M, Li Y, Glass Z, Xu Q. (2020) "Imidazole-Based Synthetic Lipidoids for In Vivo mRNA Delivery into Primary T Lymphocytes." *Angew Chem Int Ed Engl.* 59(45):20083-20089. doi: 10.1002/anie.202008082.

2) Li Y, Yang T, Yu Y, Shi N, Yang L, Glass Z, Bolinger J, Finkel IJ, Li W, Xu Q. (2018) "Combinatorial library of chalcogen-containing lipidoids for intracellular delivery of genome-editing proteins." *Biomaterials.* 178:652-662. doi: 10.1016/j.biomaterials.2018.03.011.

3) Chen J, Qiu M, Ye Z, Nyalile T, Li Y, Glass Z, Zhao X, Yang L, Chen J, Xu Q. (2021) "In situ cancer vaccination using lipidoid nanoparticles." *Sci Adv.* 7(19):eabf1244. doi: 10.1126/sciadv.abf1244.

**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 X, ZZ-ZZ-ZZ – For research use only. Not intended for diagnostic or therapeutic use.



Echelon Biosciences Inc.

Ph: 866-588-0455  
Fax: 801-588-0497

Echelon-inc.com