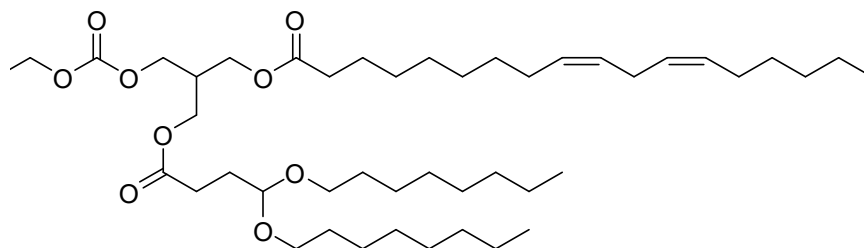


Echelon Biosciences Inc.

LP-01

Catalog number: N-1007



Molecular Formula: C₅₀H₉₃NO₉

MW: 852.29

CAS: 1799316-64-5

Alternate Name: CIN-16645, 9Z,12Z-octadecadienoic acid, 3-[4,4-bis(octyloxy)-1-oxobutoxy]-2-[[[3-(diethylamino)propoxy] carbonyl]oxy]methyl]propyl ester

Solubility: Methanol, ethanol, DMF, DMSO.

Storage and Handling: LP-01 is stable for at least 2 years when stored as an oil, protected from moisture, and light at -20°C. Solutions of LP-01 should be stored at -20°C between uses.

Background: LP-01 is an ionizable cationic lipid and has similar structure as 11-A-M (Echelon Cat#: N-1111). LP-01 can form lipid nanoparticles (LNPs) with cholesterol, DSPC, and PEG2k-DMG. The LNP nucleic acid assemblies can deliver single-guide RNA (sgRNA) targeting MHC class II transactivator (CIITA) in T cells.

References: 1) Srijani S. et al. "Compositions and methods for reducing MHC class II protein expression in a cell" WO/2022/125982. 2) Luis B. et al, "LIPIDS AND LIPID COMPOSITIONS FOR THE DELIVERY OF ACTIVE AGENTS" WO2015/95340,2015. 3) Stephen S. et al. "IONIZABLE AMINE LIPIDS AND LIPID NANOPARTICLES" WO2020/219876, 2020. 4) Lokugamage M. et al. "Constrained nanoparticles deliver siRNA and sgRNA to T cells in vivo without targeting ligands" Adv. Mater. 2019, 31(41) e1902251.

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, 2-16-24 – 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