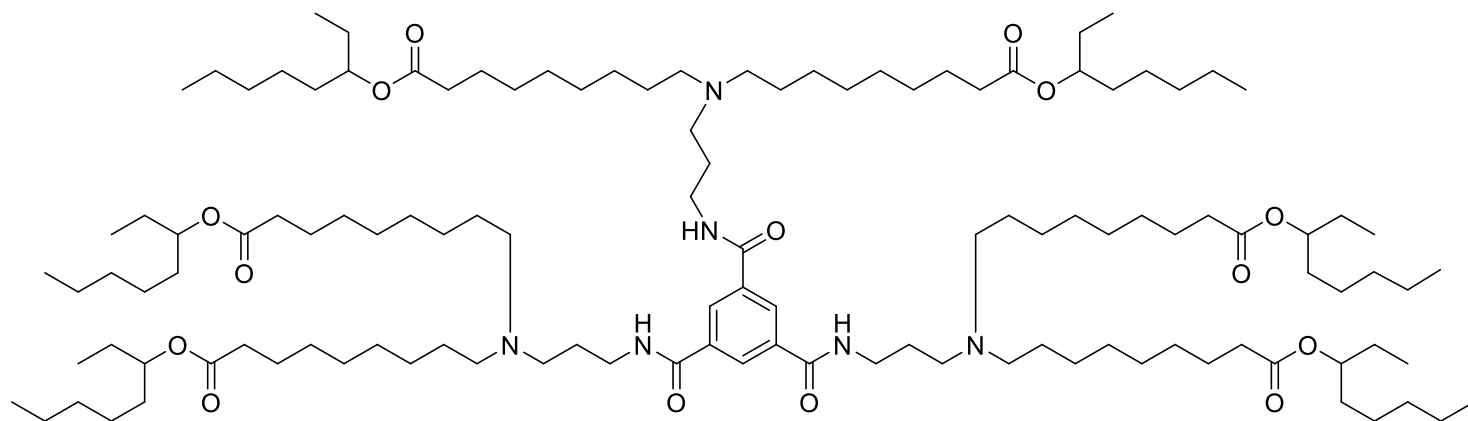


Echelon Biosciences Inc.

FTT5

Catalog number: N-1115



Molecular Formula: $C_{120}H_{222}N_6O_{15}$

MW: 1989.12

CAS: 2328129-27-5

Alternate Name: hexa(octan-3-yl) 9,9',9'',9''',9''''-((((benzene-1,3,5-tricarbonyl)tris(azanediyl))tris(propane-3,1-diyl))tris(azanetriyl))hexanonanoate

Solubility: Methanol, ethanol, DMSO.

Storage and Handling: FTT5 is stable for at least 2 years when stored as an oil, protected from moisture, and light at -20°C . Solutions of FTT5 should be stored at -20°C between uses.

Background: FTT5 (Echelon Cat#: N-1115) is a functionalized derivative of TT3 (Echelon Cat#: N-1113). FTT5 lipid like nanomaterials (LLNs) displayed two-fold potency in the delivery efficiency of mRNA to the liver, compared to TT3 LLNs. FTT5 LLNs efficient delivery long mRNAs, such as human factor VIII (hFVIII) mRNA (~4.5 kb) for expression of hFVIII protein in hemophilia A mice *in vivo*.

References: 1) Zhang X. et al. "Functionalized lipid-like nanoparticles for *in vivo* mRNA delivery and base editing". *Sci. Adv.* 2020,6(34), eabc2315. DOI: 10.1126/sciadv.abc2315

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, 3-28-22 – 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