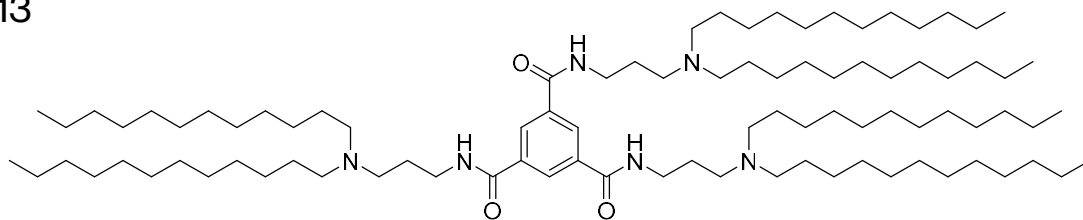


TT3

Catalog number: N-1113



Molecular Formula: C₉₀H₁₇₄N₆O₃

MW: 1388.42

CAS#: 1821214-50-9

Alternate Name: N¹,N³,N⁵-tris[3-(didodecylamino)propyl]-1,3,5-benzenetricarboxamide, O-TT3

Solubility: Methanol, ethanol, DMSO.

Storage and Handling: TT3 is stable for at least 2 years when stored as an oil, protected from moisture, and light at -20°C. Solutions of TT3 should be stored at -20°C between uses.

Background: TT3, together with helper lipid, DMG-PEG2000, Cholesterol, and mRNA formed TT3 lipid-like nanoparticles (LLNs) which improve the delivery efficiency of firefly luciferase FLuc mRNA over 350-fold *in vitro*. TT3 LLNs with human factor IX (hFIX) mRNA can efficiently improve the hFIX level to normal physiological values in FIX-knockout mice.

References: 1) Li, B. et al. "An orthogonal array optimization of lipid-like nanoparticles for mRNA delivery *in vivo*". *Nano Lett.* 2015,15(12), 8099-8107. DOI: 10.1021/acs.nanolett.5b03528

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.

