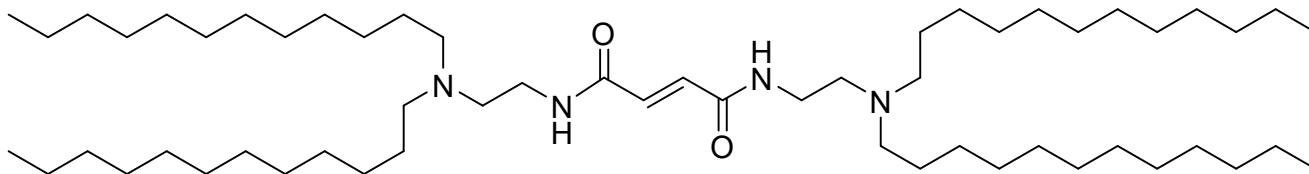


YSH-12

Catalog number: N-1009



Molecular Formula: C₅₆H₁₁₂N₄O₂

MW: 873.54

CAS#: 2959463-68-2

Alternative Names: N¹,N⁴-bis(2-(didodecylamino)ethyl)fumaramide

Solubility: chloroform, methylene chloride >20 mg/mL, ethanol 2 mg/mL with sonication

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

Background: YSH-12 is an ionizable lipid used to generate lipid nanoparticles to deliver RNA. LNPs optimized with the formula 15:10:25:2.5 (YSH-12/DOPE/cholesterol/DMG-PEG) delivered mRNA and siRNA together to Raw264.7-Luc cells. YSH-12 LNPs incorporating CRV peptide were able to target monocytes *in situ* to deliver MRSA-targeted chimeric antigen receptor (CAR) mRNA (SasA-CAR mRNA) and CASP11 siRNA. The engineered monocytes showed increased bactericidal activity in a MRSA-induced mouse model.

References: C. Tang, W. Jing, *et al.* (2024) "mRNA-Laden Lipid-Nanoparticle-Enabled *inSitu* CAR-Macrophage Engineering for the Eradication of Multidrug-Resistant Bacteria in a Sepsis Mouse Model" *ACS Nano* 18(3), 2261–2278. DOI: 10.1021/acsnano.3c10109.

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 2, 5/21/2024 – **For research use only.** Not intended for diagnostic or therapeutic use.