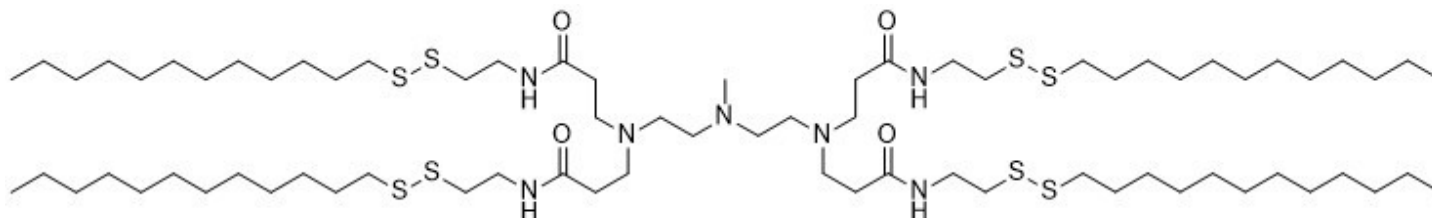


113-N16B

Catalog Number: N-1040



Molecular Formula: $C_{73}H_{147}N_7O_4S_8$

MW: 1443.5

CAS#: n/a

Solubility: methanol, ethanol, DMSO.

Storage and Handling: 113-N16B should be stored as oil at -20°C. Stock solutions of 113-N16B should be stored at -20°C between uses.

Background: 113-N16B is a biodegradable, disulfide containing ionizable lipid used in lipid nanoparticles engineered for selective mRNA delivery to the lung. 113-N16B delivered Cre mRNA primarily to lung endothelium cells after i.v. administration in a mouse model.

References: 1) Min Qiu, et al. Lung-selective mRNA delivery of synthetic lipid nanoparticles for the treatment of pulmonary lymphangioleiomyomatosis. Proc Natl Acad Sci U S A. 2022 Feb 22;119(8):e2116271119. doi: 10.1073/pnas.2116271119.

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the SDS 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.

For research use only. Not intended for diagnostic or therapeutic use.

N-1040 Technical Data Sheet / Rev 1/ 06222026