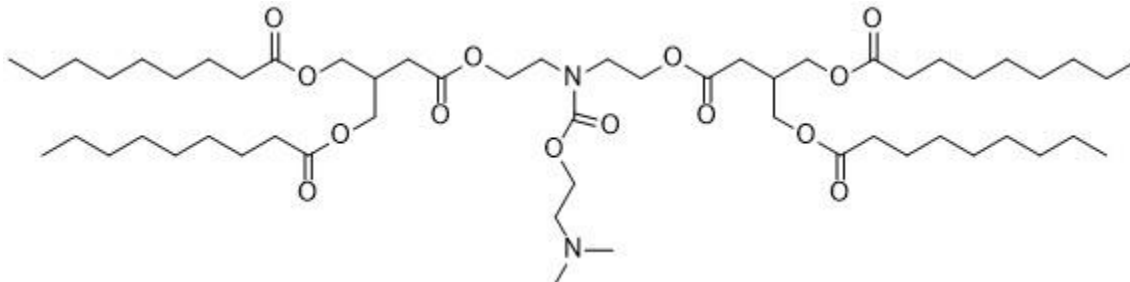


Lipid A-11 (CICL-1, L829)

Catalog Number: N-1039



Molecular Formula: C₅₅H₁₀₀N₂O₁₄

MW: 1013.41

CAS: 2996998-09-3

Alternative Names: CICL-1, L829, Lipid 829

Solubility: methanol, ethanol, DMSO.

Storage and Handling: Lipid A-11 should be stored as oil at -20°C. Stock solutions of Lipid A-11 should be stored at -20°C between uses.

Background:

To overcome the limitations of chimeric antigen receptor (CAR) T cell therapy, Capstan Therapeutics introduced a strategy using targeted lipid nanoparticles (tLNPs) to deliver mRNA to specific T cell subsets, enabling the safe and preferential generation of CAR T cells directly *in vivo*. L829-LNP showed reduced functional delivery to the liver, efficient delivery of mRNA to target immune cells, improved biodegradability, reduced reactogenicity, and improved tolerability. L829 was rapidly cleared from the liver in CD5-L829-tLNP-treated mice, whereas ALC-0315 levels remained elevated for at least 7 days.

References: 1) Theresa L. Hunter, et al. *In vivo* CAR T cell generation to treat cancer and autoimmune disease. *Science*, 2025, 388 (6753), 1311-1317

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, 05122026 – **For research use only.** Not intended for diagnostic or therapeutic use.