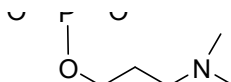


Echelon Biosciences Inc.

CAP2

Catalog number: N-1912



Molecular Formula: C₅₉H₁₀₂NO₄P

MW: 920.44

CAS: N/A

Alternate Name: bis((3S,8S,9S,10R,13R,14S,17R)-10,13-dimethyl-17-((R)-6-methylheptan-2-yl)-2,3,4,7,8,9,10,11,12,13,14,15,16,17-tetradecahydro-1H-cyclopenta[a]phenanthren-3-yl) (3-(dimethylamino)propyl) phosphate

Solubility: Ethanol, DMF, DMSO.

Storage and Handling: CAP2 is stable for at least 2 years when stored as powder, protected from moisture, and light at -20°C. Solutions of CAP2 should be stored at -20°C between uses.

Background: Lipid Nanoparticles (LNPs) made from Cholesterol-Amino-Phosphate (CAP) CAP2 can deliver traditional mRNA and self-amplifying RNA (saRNA) encoding DNA Meiotic Recombinase 1 (Dmc1) protein in spermatocytes. A single injection of CAP2 LNPs-saRNA can produce Dmc1 protein for an extended period and restore spermatid production in the Dmc1 gene knockout mouse model.

References: 1) Du S. et al. "Cholesterol-Amino-Phosphate (CAP) Derived Lipid Nanoparticles for Delivery of Self-Amplifying RNA and Restoration of Spermatogenesis in Infertile Mice" Adv. Sci. 2023, 2300188.

DOI: 10.1002/advs.202300188

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, 1-19-24 – 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