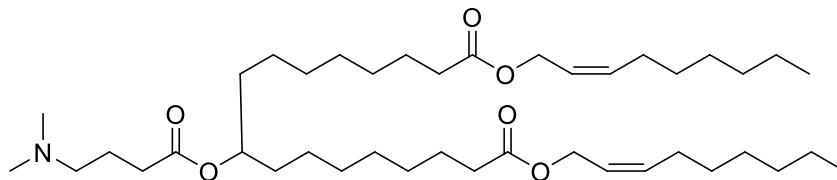


Echelon Biosciences Inc.

Lipid 319

Catalog number: N-1319



Molecular Formula: $C_{41}H_{75}NO_6$

MW: 678.05

CAS: 1351586-50-9

Alternate Name: 9-[4-(dimethylamino)-1-oxobutoxy]-heptadecanedioic acid, 1,17-di-(Z)-2-nonen-1-yl ester
Di((Z)-Non-2-en-1-yl)9-((4-(dimethylamino)butanoyl)oxy)heptadecanedioate

Solubility: Methanol, ethanol, DMSO, and dimethyl formamide.

Storage and Handling: Lipid 319 is stable for at least one year when stored as a solid, protected from moisture, and light at -20°C . Solutions of Lipid 319 should be stored at -20°C between uses.

Background: Lipid 319 is an ionizable and biodegradable lipid. Lipid 319 nanoparticle has better delivery efficacy and biodegradability than MC3 in the delivery of mRNA in vivo. MC3 is the key delivery nanoparticle component of the first siRNA drug Onpattro.

References: 1) Maier M. et al. "Biodegradable Lipids Enabling Rapidly Eliminated Lipid Nanoparticles for Systemic Delivery of RNAi Therapeutics." *Mol Ther.* 2013, *21(8)*,1570-1578
<https://www.ncbi.nlm.nih.gov/pmc/articles/PMC3734658/>

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, 3-28-22 – 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