

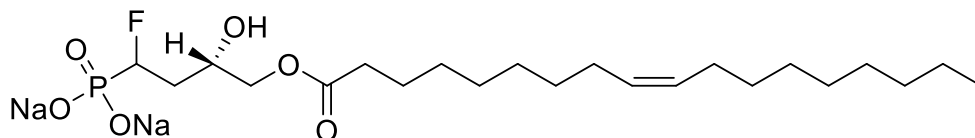
(3S) 1-Fluoro-3-hydroxy-4-(oleoyloxy)butyl-1-phosphonate (XY-17)

Catalog number: L-9118

Molecular Formula: $C_{22}H_{40}FNa_2O_6P$

MW: 790.50

CAS: 565438-23-5



Solubility: Methanol, > 1 mg/mL. Stock solutions can be further diluted into buffer.

Storage and Handling: Store dry at $-20\text{ }^{\circ}\text{C}$. Stock solutions should be stored frozen ($-20\text{ }^{\circ}\text{C}$ or below). Storage in basic buffers (pH > 9.0) and acidic buffers (pH < 4.0) will result in slow decomposition.

Background: XY-17 is a fluorinated phosphonate analog of lysophosphatidic acid (LPA). XY-17 is not hydrolyzed by lipid phosphate phosphatase (LPP) and acts as an agonist of the LPA3 receptor.

References: 1) Xu, Y., M. Tanaka, et al. (2004). "Alkyl lysophosphatidic acid and fluoromethylene phosphonate analogs as metabolically-stabilized agonists for LPA receptors." *Bioorg Med Chem Lett* 14(21): 5323-8.

2) Zhao, Y., P. V. Usatyuk, et al. (2005). "Lipid phosphate phosphatase-1 regulates lysophosphatidic acid-induced calcium release, NF-kappaB activation and interleukin-8 secretion in human bronchial epithelial cells." *Biochem. J.* 385(2): 493-502.

3) Prestwich, G., Y. Xu, et al. (2005). "New metabolically stabilized analogues of lysophosphatidic acid: agonists, antagonists and enzyme inhibitors." *Biochem Soc Trans* 33: 1357-1361.

4) W.N. Ayre, T. Scott, et al. (2016) "Fluorophosphonate-functionalised titanium via a pre-adsorbed alkane phosphonic acid: a novel dual action surface finish for bone regenerative applications." *J Mater Sci Mater Med.* Feb;27(2):36. doi: 10.1007/s10856-015-5644-7

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 4, 2-11-19 – 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