



Echelon Biosciences Inc.
675 Arapleen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

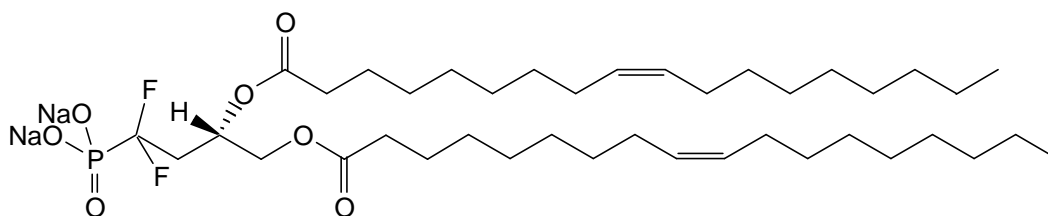
Technical Data Sheet

For research use only

*Not intended or approved for
diagnostic or therapeutic use.*

Product Name: (3S)-[1,1-Difluoro-3,4-bis(oleoyloxy)butyl]phosphonate (Sodium Salt) [1,2-Dioleoyl-3-CF₂-PA, XY-14]

Product Number: L-9218



Formula: C₄₀H₇₁F₂Na₂O₇P

MW : 778.94

CAS: 478529-66-7

Solubility: 2:1 DMSO:H₂O, 1 mg/mL

Specifications and Storage Recommendations: Phospholipids/lysophospholipids are fairly stable at room temperature in dry form. They are best stored dry at -80°C. Plastic containers may be used for the storage of short-chain phospholipids; however, longer-chain phospholipids should be stored in glass containers to prevent material loss due to adsorption on the vessel surface. Storage in basic buffers (pH > 9.0) and acidic buffers (pH < 4.0) will result in slow hydrolysis and migration of the ester chains and may cause other forms of decomposition.

Field of Interest: (3S)-[1,1-Difluoro-3,4-bis(oleoyloxy)butyl] phosphonate is a potent competitive inhibitor of Lipid phosphate phosphatase 1 (LPP1), inhibits the dephosphorylation of LPA by intact platelets, and potentiates platelet aggregation and shape change responses to LPA.

Reference: S.S. Smyth *et al.* "Lipid Phosphate Phosphatases Regulate Lysophosphatidic Acid Production and Signaling in Platelets" *J. Biol. Chem.* **2003**, 278, 43214-43223.

Hazardous Properties and Cautions: No hazardous or toxic properties are known for this substance. (3S)-[1,1-Difluoro-3,4-bis(oleoyloxy)butyl] phosphonate (Sodium Salt) is not listed as a hazardous waste or as a toxic chemical subject to release reporting under the Emergency Planning and Community Right-to-Know Act. For solutions containing methanol or other solvents, see MSDS for phospholipids in solution, available upon request. (3S)-[1,1-Difluoro-3,4-bis(oleoyloxy)butyl] phosphonate (Sodium Salt) is not listed in the Chemical Inventory of the Toxic Substance Control Act, and 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.