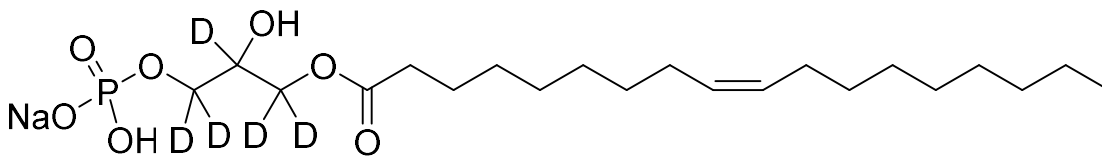


Echelon Biosciences Inc.

1-O-Oleoyl-glycerol-3-phosphate-d5 (Oleoyl LPA d5)

Catalog number: L-0181d



Molecular Formula: C₂₁H₃₅D₅NaO₇P

MW: 463.54

CAS: n/a

Alternative Names: n/a

Solubility: ~ 5mg/mL in H₂O

Storage and Handling: 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

Background: Lysophosphatidic Acid (LPA) is a small lysophospholipid involved in diverse cellular processes such as cell proliferation, chemotaxis, platelet aggregation, wound healing, angiogenesis, tumor invasion, and smooth muscle contraction. LPA binds to several G-coupled protein receptors to initiate its biological functions. In cancer, LPA primarily promotes cell survival, migration and invasion.

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, 09-23-21 – 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