

Echelon Biosciences Inc.

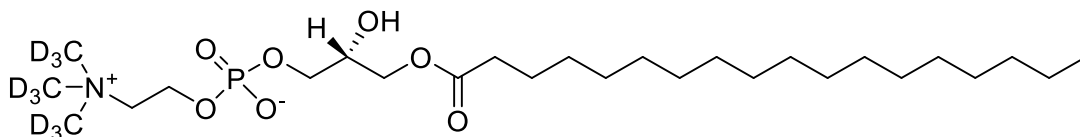
S-LysoPC-d₉ (1-Stearoyl-2-hydroxy-sn-glycero-3-phosphocholine-d₉)

Catalog number: L-1518d

Molecular Formula: C₂₆H₄₅D₉NO₇P

MW: 532.75

CAS: n/a



Alternative Names: 18:0 LPC

Solubility: chloroform, methanol

Storage and Handling: Store dry at -20 °C. Stock solutions should be stored frozen (-20 °C or below)

Background: LysoPC is generated through hydrolysis of the sn-2 acyl chain of PC by phospholipase A2. It is a major component of oxidized low-density lipoprotein and is believed to have an important role in inflammatory diseases and atherosclerosis. LysoPC is the substrate for Autotaxin/LysoPLD generating lysophosphatidic acid which has an important role in cell motility. This analog is deuterated on the choline headgroup.

References: Matsumoto, T., Kobayashi, T., Kamata, K. (2007) "Role of lysophosphatidylcholine (LPC) in atherosclerosis", *Curr. Med. Chem.* 14, 3209-3220.

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, 5-12-20 – 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