

(*R,R*)-2,2'-Bisdodecyl-LBPA (C₁₂-ether LBPA)

Catalog number: L-B120e

Molecular Formula: C₃₀H₆₆NO₈P

MW: 599.82

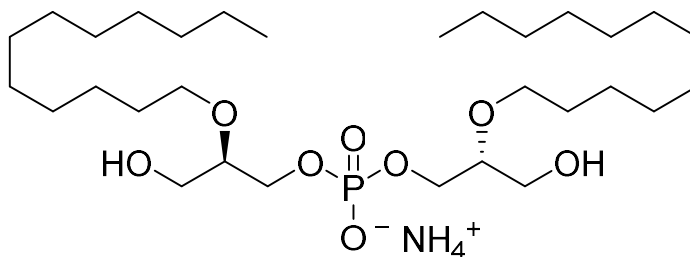
CAS#: n/a

Solubility: MeOH:CHCl₃ (1:1 v/v) ~ 1mg/mL; *t*-BuOH: water (6:1 v/v) ~ 1mg/mL

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

Background: Lysobisphosphatidic acids (LBPA)s, known also as bis-(monoacylglycerol)phosphates (BMPs), are specialized lipids reported to play a role in intracellular protein and lipid transport in healthy cells. Their accumulation was observed in pathologic liver tissue, and they also serve as antigens for auto-antibody generation in a human autoimmune condition termed antiphospholipid syndrome. Accumulation of LBPA in intracellular, often multilamellar membranes is related to biomembrane polymorphism which may impact intracellular cholesterol transport. C12-Ether LBPA is an ether analog of the naturally occurring, biologically active isomer of LPBA.

References: G. Jiang, *et al.* (2005) "Concise Synthesis of Ether Analogues of Lysobisphosphatidic Acid" *Org. Lett.* 7(18): 3837-3840.



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 2, 6/27/2024 – **For research use only.** Not intended for diagnostic or therapeutic use.