



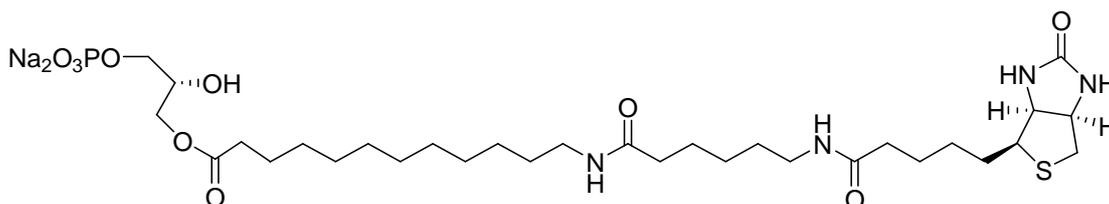
Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only
*Not intended or approved for
diagnostic or therapeutic use.*

Product Name: LPA-biotin

Product Number: L-012B



Formula: C₃₁H₅₇N₄O₁₀PS, **MW:** 708.85

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: Many bioactive lysophospholipid compounds exhibit pleiomorphic effects on numerous cell lineages including ovarian cancer cells as described in many publications. Depending on the application, choices can be made from a variety of lysophospholipid analogs including saturated or polyunsaturated lipid species, non-migrating and non-hydrolyzable analogs, labeled derivatives, and metabolically stable agonists and antagonists.

Hazardous Properties and Cautions: No hazardous or toxic properties are known for this substance. LPA-biotin 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. LPA-biotin 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.

Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications and methods and use, may be covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our trademarks and technologies at busdev@echelon-inc.com.

TDS L-012B Rev: 1 (01/24/05)