

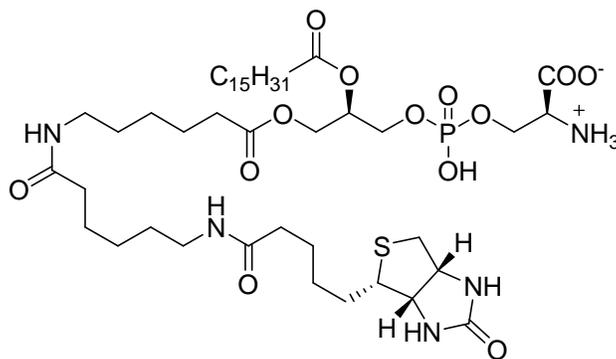


Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only

Product Name: **Biotin-Labeled Phosphatidylserine**



Catalog Number	MW (g/mol)	Chemical Formula	Solubility
L-31B16	950.17	C ₄₄ H ₈₀ N ₅ O ₁₃ PS	MeOH, EtOH, CHCl ₃ >1 mg/ml

Storage: Biotin-Labeled Phosphatidylserine (biotin-PS) is stable for at least one year when stored as a solid, protected from moisture, and light at -20 °C. Reconstitute with methanol and dilute with water or neutral pH, buffered salt solutions, i.e. PBS, TBS, etc. Should be stored in glass containers to prevent material loss due to adsorption to the vessel surface. Storage in basic (pH > 9) or acidic (pH < 4) buffers will result in slow decomposition of the product. After reconstitution, solutions of biotin-PS should be flash frozen in liquid nitrogen and stored at -20 °C between uses. Biotin-PS is stable for at least three months when handled in this way. Repeated freeze/thaw cycles do not affect biotin-PS. Do not store reconstituted biotin-PS, at 4°C for more than 2-3 days.

Field of Interest: Phosphatidylserine (PS) is a negatively charged phospholipid located in the inner leaflet of the cell membrane. During apoptosis PS is no longer restricted to the inner leaflet of the plasma membrane and migrates to the outer leaflet. Recognition of extra-leaflet PS by Anxin V is the basis of popular apoptosis assays. Similarly, during platelet activation, PS is translocated to the outer side where it facilitates assembly of the prothrombinase complex.

Hazardous Properties and Cautions: No hazardous or toxic properties are known for this substance. Biotin-Labeled Phosphatidylserine 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 Phosphoinositides in Solution available on request. Biotin-Labeled Phosphatidylserine is not listed on 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.

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.

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.