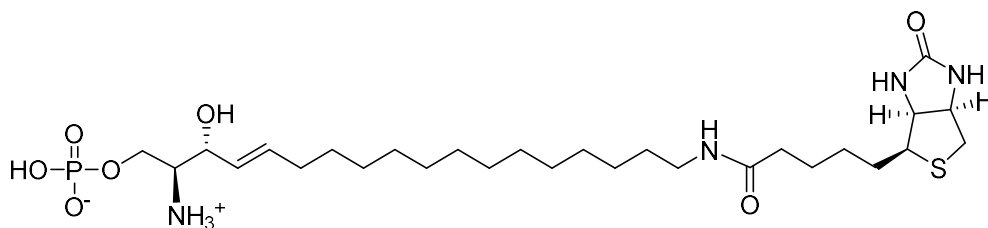


Sphingosine 1-phosphate biotin (S1P-biotin)

Catalog number: S-200B



Molecular Formula: C₂₈H₅₃N₄O₇PS

MW: 620.8

CAS#: 799812-63-8

Solubility: 1:1 H₂O:DMSO, 1 mg/ml

Storage and Handling: Sphingosine 1-phosphate biotin is relatively stable at room temperature as a solid. It is best stored as a dried solid, protected from light at -20°C. Solutions/emulsions of sphingosines should be kept at -20°C. Storage in basic buffers (pH > 9.0) or acidic buffers (pH < 4.0) may cause decomposition.

Background: Sphingosine 1-phosphate (S1P) is a bioactive lysophospholipid formed by phosphorylation of sphingosine by sphingosine kinases. S1P binds to five G-protein coupled receptors designated S1P₁₋₅ leading to cell proliferation, motility, and has a role in regulation cell shape. In addition, S1P acts as an intracellular mediator in a receptor-independent manner regulating cell survival and Ca²⁺ homeostasis. This S1P analog is labeled with biotin at the 3-position and can be bound to streptavidin coated surfaces such as beads, microtiter plates, etc.

References: 1) N. Ieronimakis, *et al.* (2013) "Increased sphingosine-1-phosphate improves muscle regeneration in acutely injured mdx mice" *Skelet. Muscle* 3(1): 20.

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 3, 5/13/24– **For research use only.** Not intended for diagnostic or therapeutic use.