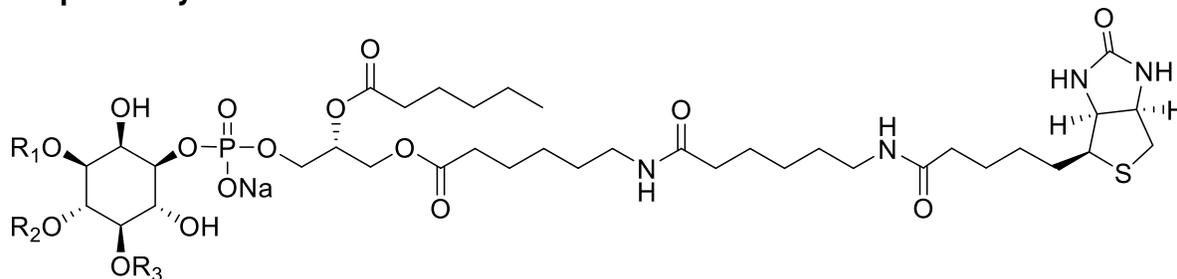


Echelon Biosciences Inc.

Biotin Phosphatidylinositols



| Catalog Number | R ₁ | R ₂ | R ₃ | MW (g/mol) | Chemical Formula |
|----------------|---------------------------------|---------------------------------|---------------------------------|------------|---|
| C-00B6 | H | H | H | 907.0 | C ₃₇ H ₆₄ N ₄ NaO ₁₆ PS |
| C-03B6 | PO ₃ Na ₂ | H | H | 1030.9 | C ₃₇ H ₆₃ N ₄ Na ₃ O ₁₉ P ₂ S |
| C-04B6 | H | PO ₃ Na ₂ | H | 1030.9 | C ₃₇ H ₆₃ N ₄ Na ₃ O ₁₉ P ₂ S |
| C-05B6 | H | H | PO ₃ Na ₂ | 1030.9 | C ₃₇ H ₆₃ N ₄ Na ₃ O ₁₉ P ₂ S |
| C-34B6 | PO ₃ Na ₂ | PO ₃ Na ₂ | H | 1154.8 | C ₃₇ H ₆₂ N ₄ Na ₅ O ₂₂ P ₃ S |
| C-35B6 | PO ₃ Na ₂ | H | PO ₃ Na ₂ | 1154.8 | C ₃₇ H ₆₂ N ₄ Na ₅ O ₂₂ P ₃ S |
| C-45B6 | H | PO ₃ Na ₂ | PO ₃ Na ₂ | 1154.8 | C ₃₇ H ₆₂ N ₄ Na ₅ O ₂₂ P ₃ S |
| C-39B6 | PO ₃ Na ₂ | PO ₃ Na ₂ | PO ₃ Na ₂ | 1278.8 | C ₃₇ H ₆₁ N ₄ Na ₇ O ₂₅ P ₄ S |

Solubility: 5mg/mL in H₂O

Storage and Handling: Phosphatidylinositol polyphosphates (PtdInsP_ns) and analogs are stable for at least one year when stored as a solid, protected from moisture, at -20 °C or below. Storage in basic (pH > 9) or acidic (pH < 4) buffers may cause decomposition. After reconstitution, solutions of PtdInsP_ns should be stored at -20 °C or below. PtdInsP_ns are stable for at least three months when handled in this way. Repeated freeze/thaw cycles do not affect PtdInsP_ns. Do not store reconstituted PtdInsP_ns at 4 °C for more than 2-3 days.

Background: These water-soluble phosphatidylinositide analogs are labeled with biotin at the sn-1 position and can be bound to streptavidin coated surfaces such as beads, microtiter plates, etc.

References: C-00B6: Bidlingmaier, S., Y. Wang, et al. (2011). "Comprehensive analysis of yeast surface displayed cDNA library selection outputs by exon microarray to identify novel protein-ligand interactions." *Mol Cell Proteomics* 10(3): M110 005116. C-03B6: Jank, T., K. E. Böhmer, et al. (2012). "Domain organization of Legionella effector SetA." *Cellular Microbiology* 14(6): 852. C-39B6: Anzelon, A. N., H. Wu, et al. (2003). "Pten inactivation alters peripheral B lymphocyte fate and reconstitutes CD19 function." *Nat Immunol* 4(3): 287-94. C-45B6: A. L. Egea-Jimenez, R. Gallardo, A. Garcia-Pino, Y. Ivarsson, A. M. Wawrzyniak, R. Kashyap, R. Loris, J. Schymkowitz, F. Rousseau and P. Zimmermann. (2016) "Frizzled 7 and PIP2 binding by syntenin PDZ2 domain supports Frizzled 7 trafficking and signalling" *Nat Commun* 7, DOI: 10.1038/ncomms12101. *See the product webpages for additional references.*

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 6, 10-31-18 – 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