



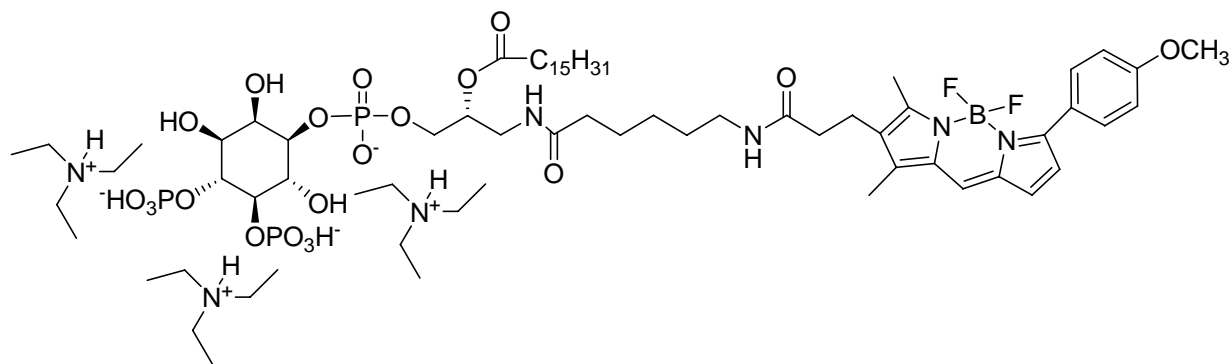
Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108 Telephone
866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only

Not intended or approved for
diagnostic or therapeutic use.

Product Name **Bodipy®-TMR-PI(4,5)P₂, C₁₆**



Product Number: C-45M16a **Formula:** C₇₀H₁₂₇BF₂N₇O₂₀P₃ (triethylammonium salt) **MW:** 1528.5

Solubility: >5mg/mL in H₂O, <5 mg/mL CHCl₃

Absorbtion/emission: 544/574 nm

Specifications and Storage Recommendations: Protect from light. Phosphoinositide polyphosphates (PtdInsP_ns) are relatively stable at room temperature as solids. Solutions of short-chain and labeled PtdInsP_ns should be stored in plastic containers at 4°C or may be flash frozen in liquid nitrogen for prolonged storage at -80°C. Storage in basic buffers (pH > 9.0) will result in slow hydrolysis of the ester chain, and may cause phosphate migration to occur. Storage in acidic buffers (pH < 4.0) may cause decomposition or phosphate migration.

Field Of Interest: PtdInsP_ns compounds are employed as substrates for kinases, phosphatases and binding proteins as described in many publications. Depending on the application, the lyophilized powders may be dissolved or suspended in water, buffers, or mixed solvents such as CHCl₃-MeOH and others employed for preparing mixed phospholipid liposomes.

Hazardous Properties and Cautions: No hazardous or toxic properties are known for this substance. BODIPY TMR-PtdInsP_ns are 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. BODIPY TMR-PtdInsP_ns are 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.