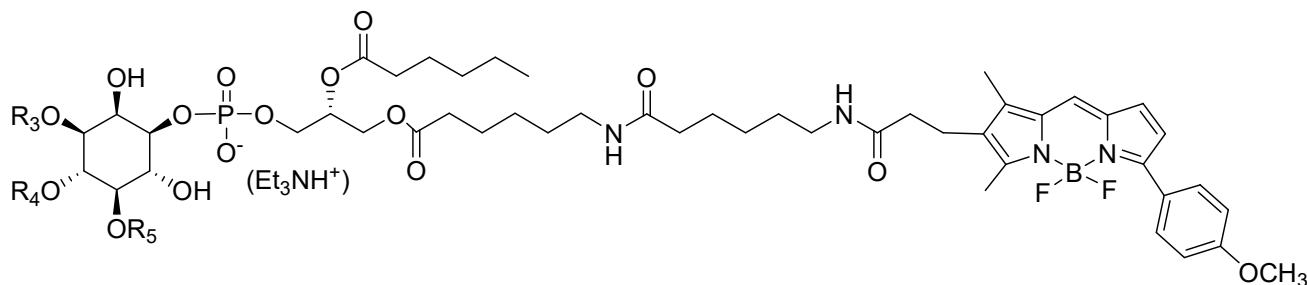


BODIPY® TMR Phosphatidylinositols



Catalog Number	R ₃	R ₄	R ₅	MW (g/mol)	Chemical Formula
C-00M6	H	H	H	1140.06	C ₅₄ H ₈₅ BF ₂ N ₅ O ₁₆ P
C-03M6	PO ₃ (HNEt ₃) ₂	H	H	1422.42	C ₆₆ H ₁₁₆ BF ₂ N ₇ O ₁₉ P ₂
C-04M6	H	PO ₃ (HNEt ₃) ₂	H	1422.42	C ₆₆ H ₁₁₆ BF ₂ N ₇ O ₁₉ P ₂
C-05M6	H	H	PO ₃ (HNEt ₃) ₂	1422.42	C ₆₆ H ₁₁₆ BF ₂ N ₇ O ₁₉ P ₂
C-34M6	PO ₃ (HNEt ₃) ₂	PO ₃ (HNEt ₃) ₂	H	1704.78	C ₇₈ H ₁₄₇ BF ₂ N ₉ O ₂₂ P ₃
C-35M6	PO ₃ (HNEt ₃) ₂	H	PO ₃ (HNEt ₃) ₂	1704.78	C ₇₈ H ₁₄₇ BF ₂ N ₉ O ₂₂ P ₃
C-45M6	H	PO ₃ (HNEt ₃) ₂	PO ₃ (HNEt ₃) ₂	1704.78	C ₇₈ H ₁₄₇ BF ₂ N ₉ O ₂₂ P ₃
C-39M6	PO ₃ (HNEt ₃) ₂	PO ₃ (HNEt ₃) ₂	PO ₃ (HNEt ₃) ₂	1987.14	C ₉₀ H ₁₇₈ BF ₂ N ₁₁ O ₂₅ P ₄

Solubility: 5mg/mL in H₂O

Excitation/Emission: 542/574 nm

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. Longer-chain PtdInsP_ns should be stored in glass containers to prevent material loss due to absorption to the vessel surface. Storage in basic (pH > 9) or acidic (pH < 4) buffers may cause decomposition. Brief sonication and heating (~50 °C) can help longer chain PtdInsP_ns solubilize. 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 the fluorophore Bodipy® TMR at the sn-1 position.

References: **C-00M6** Stankewicz, C.& Rininsland, F. H. (2016) A robust screen for inhibitors and enhancers of phosphoinositide-3 kinase (PI3K) activities by ratiometric fluorescence superquenching. *J. Biomol. Screen.*11(4), 413-422. **C-45M6:** Cabanos C, Wang M, Han X, Hansen SB (2017) A Soluble Fluorescent Binding Assay Reveals PIP2 Antagonism of TREK-1 Channels. *Cell Rep* 20(6):1287–1294, See 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 7, 5/2/24– **For research use only.** Not intended for diagnostic or therapeutic use.