

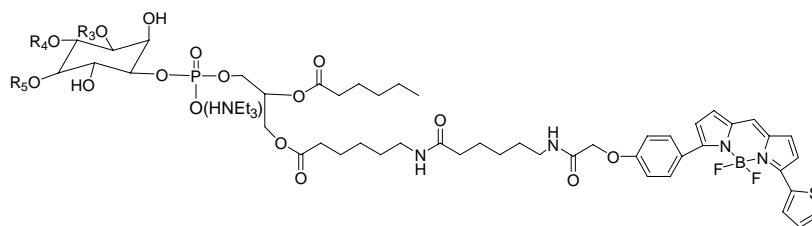


Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108 Telephone
866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only
Not intended or approved for
diagnostic or therapeutic use.

Product Name: BODIPY[®] TR Probes (589/617), C₆



Structure of BODIPY[®] TR-PtdIns; R = H or PO₃(HNEt₃)₂

Product	R ₃	R ₄	R ₅	Chemical Formula	MW (g/mol)
C-00R6	H	H	H	C ₅₄ H ₇₉ BF ₂ N ₅ O ₁₆ PS	1166.08
C-03R6	PO ₃ (HNEt ₃) ₂	H	H	C ₆₆ H ₁₁₀ BF ₂ N ₇ O ₁₉ P ₂ S	1448.44
C-04R6	H	PO ₃ (HNEt ₃) ₂	H	C ₆₆ H ₁₁₀ BF ₂ N ₇ O ₁₉ P ₂ S	1448.44
C-05R6	H	H	PO ₃ (HNEt ₃) ₂	C ₆₆ H ₁₁₀ BF ₂ N ₇ O ₁₉ P ₂ S	1448.44
C-34R6	PO ₃ (HNEt ₃) ₂	PO ₃ (HNEt ₃) ₂	H	C ₇₈ H ₁₄₁ BF ₂ N ₉ O ₂₂ P ₃ S	1730.80
C-35R6	PO ₃ (HNEt ₃) ₂	H	PO ₃ (HNEt ₃) ₂	C ₇₈ H ₁₄₁ BF ₂ N ₉ O ₂₂ P ₃ S	1730.80
C-45R6	H	PO ₃ (HNEt ₃) ₂	PO ₃ (HNEt ₃) ₂	C ₇₈ H ₁₄₁ BF ₂ N ₉ O ₂₂ P ₃ S	1730.80
C-39R6	PO ₃ (HNEt ₃) ₂	PO ₃ (HNEt ₃) ₂	PO ₃ (HNEt ₃) ₂	C ₉₀ H ₁₇₂ BF ₂ N ₁₁ O ₂₅ P ₄ S	2013.16

SOLUBILITY: < 5 mg/mL in H₂O

STORAGE: Phosphatidylinositol polyphosphates (PtdInsP_ns) and analogs are stable for at least one year when stored as a solid, protected from moisture, at -20 °C. Reconstitute with water or neutral pH, buffered salt solutions, i.e. PBS, TBS, etc. Storage in basic buffers (pH > 9) will result in slow hydrolysis of the ester chains, and may cause phosphate or acyl migration to occur. Storage in acidic buffers (pH < 4) may cause decomposition or phosphate migration. After reconstitution, solutions of PtdInsP_ns should be flash frozen in liquid nitrogen and stored at -20 °C between uses. 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.

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

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.

AND CAUTIONS:

No hazardous or toxic properties are known for this substance. C6-BODIPY TR-PtdInsP_ns 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. C6-BODIPY TR-PtdInsP_ns 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.