

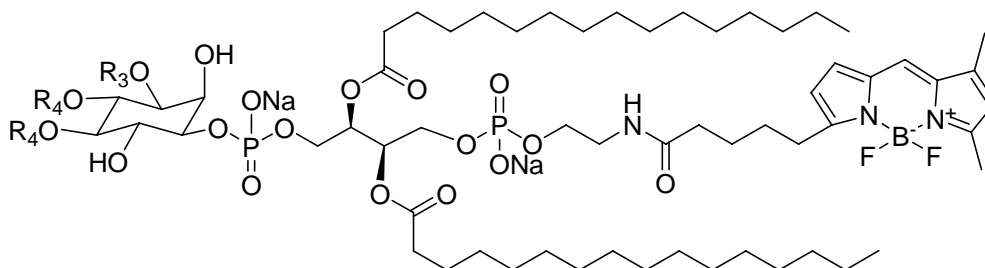


Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108 Telephone
866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

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

Product Name: Bodipy-FL Hybrid-Phosphoinositides



Product	R ₃	R ₄	R ₅	Chemical Formula	MW (g/mol)
H-34FL	PO ₃ Na ₂	PO ₃ Na ₂	H	C ₆₀ H ₁₀₀ BF ₂ N ₃ Na ₆ O ₂₄ P ₄	1558.1
H-45FL	H	PO ₃ Na ₂	PO ₃ Na ₂	C ₆₀ H ₁₀₀ BF ₂ N ₃ Na ₆ O ₂₄ P ₄	1558.1
H-39FL	PO ₃ Na ₂	PO ₃ Na ₂	PO ₃ Na ₂	C ₆₀ H ₉₉ BF ₂ N ₃ Na ₈ O ₂₇ P ₅	1682.0

Storage:

Hybrid Phosphatidylinositol polyphosphates (PtdInsP_ns) and analogs are stable for at least one year when stored as a solid, protected from moisture, at -20 °C. Hybrid PtdInsP_ns should be stored in glass containers or low-binding polypropylene tubes to prevent material loss due to absorption to the vessel surface. Storage in basic solutions (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.

Reconstitution:

Reconstitute with water or neutral pH, buffered salt solutions, i.e. PBS, TBS, etc. Hybrid PtdInsP_ns have very limited solubility in CHCl₃-MeOH and other organic solvents, and are not recommended for preparing phospholipid liposomes.

Suggested Use:

Phosphoinositides and synthetic analogs are employed as substrates for kinases, phosphatases, and binding proteins as

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.

described in many publications. Other *in vitro* and cellular applications are possible, but have not been verified by Echelon Biosciences. Please check our web site, www.echelon-inc.com, for updated technical or product application information; or call our customer service department at 1-866-588-0455.

Reference:

P.W. Rzepecki and G.D. Prestwich (2002) Synthesis of hybrid lipid probes: derivatives of phosphatidylethanolamine-extended phosphatidylinositol 4,5-bisphosphate (Pea-PIP₂). *J Org Chem*, **67**, 5454-60.