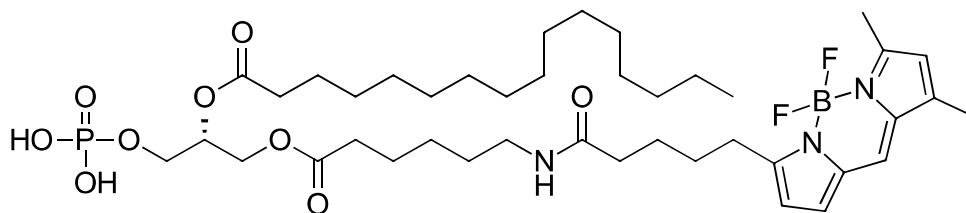


Phosphatidic Acid-Bodipy-FL (PA-Bodipy-FL)

Catalog Number: L-41F16



Molecular Formula: C₄₁H₆₇BF₂N₃O₉P

MW: 825.77

CAS#: n/a

Alternative Names: Phosphatidic Acid-Bodipy-FL, 6:0/16:0 PA-Bodipy-FL

Solubility: >5 mg/mL CHCl₃, CH₂Cl₂, 1 mg/mL MeOH

Excitation/Emission: 505/520 nm

Storage and Handling: Store dry at -20 °C protected from light. Stock solutions should be stored frozen (-20 °C or below).

Background: Phosphatidic acid (PA) is a minor component of cell membranes, constituting 1 – 2% of all phospholipids in humans. PA is made by phospholipase D from phosphatidylcholine and is used as a substrate for synthesis of other phospholipids. PA is implied to play a role in both vesicle fusion and fission. This analog is labeled on the *sn*-1 acyl chain with the green fluorophore Bodipy-FL.

References: 1) Wang, X., Devaiah, S.P., Zhang, W., Welte, R. "Signaling functions of Phosphatidic Acid" (2006), Prog. Lipid Res. 45: 250-278.
2) Foster, D.A., et al. (2014) "Phospholipase D and the Maintenance of Phosphatidic Acid Levels for Regulation of mTOR" J. Biol. Chem. 289: 22583-22588.
3) Raben, D.M., Barber, C.N. (2017) "Phosphatidic acid and neurotransmission" Adv. Biol. Regul. 63:15-21.

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the SDS 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.

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

L-41F16 Technical Data Sheet / Rev #1 / 6/10/2026