

FS-3 (LysoPLD/ATX substrate)

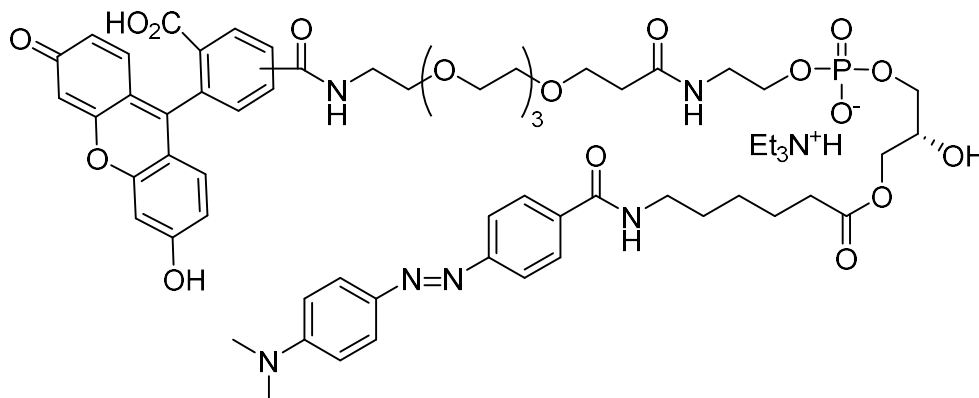
Catalog number: L-2000

Molecular Formula: C₆₄H₈₄N₇O₁₉P

MW: 1286.4

CAS: 889132-42-7

Solubility: 5 mg/mL in H₂O or MeOH



Storage and Handling: Store dry at -20 °C protected from light. Stock solutions should be stored frozen (-20 °C or below). Reconstitute with water or neutral pH, buffered salt solutions, i.e. PBS, TBS, etc. Storage in basic (pH > 9) or acidic (pH < 4) buffers may cause decomposition. Do not store reconstituted at 4°C for more than 2-3 days

Background: Autotaxin which has lysophospholipase D (lysoPLD) activity, cleaves choline from lysophosphatidylcholine forming lysophosphatidic acid (LPA), a potent mitogen that has been implicated in the pathophysiology of ovarian cancer. LysoPLD/ATX has been demonstrated to increase cell motility, neovascularization, proliferation, and aggressiveness of tumors and is upregulated in numerous cancer lineages (non-small cell lung, glioma, mammary carcinoma, renal cell carcinoma, hepatocellular carcinoma). In addition, dysregulation of the ATX/LPA pathway is central to the pathophysiology of idiopathic pulmonary fibrosis, rheumatoid arthritis, and other inflammatory diseases. Modulation of ATX activity through small-molecules is a currently underexplored target, but one with high potential for novel cancer and anti-inflammatory therapeutics.

Application: The substrate can be used in the typical conditions for *in vitro* Autotaxin assays (e.g. 50 mM Tris, 140 mM NaCl, 5 mM KCl, 1 mM CaCl₂, 1 mM MgCl₂, pH 8.0). Substrate concentrations can vary from 0.5-5 μM and incubation times (37°C) can vary from a few hours to overnight depending on enzyme activity and concentration. When using the substrate with biological samples (e.g. plasma, serum, ascites...), it is advisable to dilute said sample 2-10X in PBS. The excitation and emission spectra of FS-3 are typical of fluorescein, with maxima at about 494 nm and 520 nm respectively. At higher substrate concentrations, it may be necessary to use shorter excitation and longer emission wavelengths.

References: 1) Ferguson, C. G., C. S. Bigman, et al. (2006). "Fluorogenic phospholipid substrate to detect lysophospholipase D/autotaxin activity." *Org Lett* 8(10): 2023-6.

2) Hoeglund, A. B., H. E. Bostic, et al. (2010). "Optimization of a pipemidic acid autotaxin inhibitor." *J Med Chem* 53(3): 1056-66.

See website 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.