

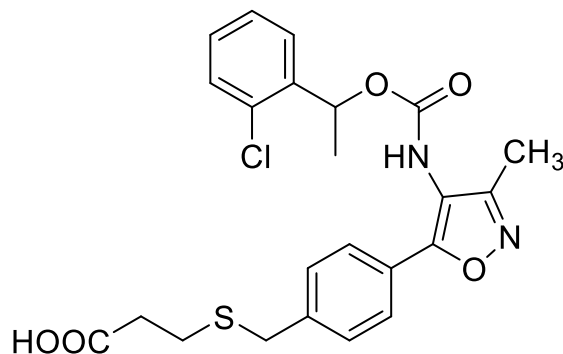
Ki16425 (LPA_{1/3} Antagonist)

Catalog number: B-0730

Molecular Formula: C₂₃H₂₃ClN₂O₅S

MW: 474.10

CAS: 355025-24-0



Alternate Name: 3-[[[4-[4-[[[1-(2-Chlorophenyl)ethoxy]carbonyl]amino]-3-methyl-5-isoxazolyl]phenyl]methyl]-thio]propanoic acid

Solubility: DMSO, > 5 mg/mL

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

Background: To date, six types of receptors, LPA₁₋₆ have been identified and are responsible for most of the biological activities of LPA. Recent studies on gene targeting in mice and family diseases of these receptors revealed that LPA is involved in various patho-physiological states. LPA antagonists have attracted considerable attention and numerous small molecules having LPA antagonistic activity have been reported. Ki16425 is an antagonist of LPA₁ and LPA₃ receptors with moderate activity against LPA₂.

References: 1). Hideo Ohta, et al “Ki16425, a Subtype-Selective Antagonist for EDG-Family Lysophosphatidic Acid Receptors” *Molecular Pharmacology*, 2003, 166(4), 994-1005; 2). Jing Zhao, et al “Lysophosphatidic Acid Receptor 1 antagonist Ki16425 Blunts Abdominal and Systemic Inflammation in a Mouse Model of Peritoneal Sepsis” *Translational Research*, 2015, 166(1), 80-88

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.

