

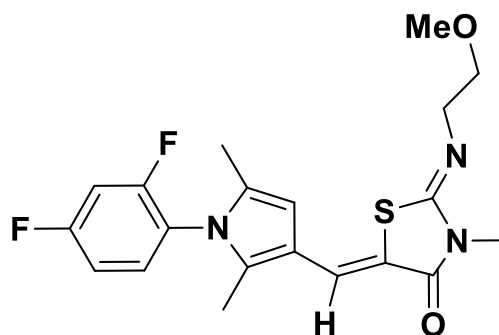
CYM50308 (S1P₄ Agonist)

Catalog number: B-0060

Molecular Formula: C₂₀H₂₁F₂N₃O₂S

MW: 405.46

CAS: 1345858-76-5



Alternate Names: (2Z,5Z)-5-[[1-(2,4-Difluorophenyl)-2,5-dimethyl-1H-pyrrol-3-yl]methylene]-2-[(2-methoxyethyl)imino]-3-methyl-4-thiazolidinone, ML248

Solubility: up to 10 mM in DMSO

Storage and Handling: Store dry at 4 °C. Stock solutions should be stored frozen (-20 °C or below).

Background: CYM50308 is a potent and selective S1P₄ receptor agonist with EC₅₀ of 37.7 - 79.1 nM. It is inactive against the S1P₁, S1P₂, and S1P₃ receptors up to 25 μM and the S1P₅ receptor up to 2.1 μM.

References: 1) Urbano *et al.* Discovery, synthesis and SAR analysis of novel selective small molecule S1P₄-R agonists based on a (2Z,5Z)-5-((pyrrol-3-yl)methylene)-3-alkyl-2-(alkylimino)thiazolidin-4-one chemotype, *Bioorganic & Medicinal Chemistry Letters*, 2011, 6739-6745. 2) Guerrero *et al.* Identification of a novel agonist of the sphingosine 1-phosphate receptor 4 (S1P₄), Probe Reports from the NIH Molecular Libraries Program. <http://www.ncbi.nlm.nih.gov/books/NBK143556/>

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.

