

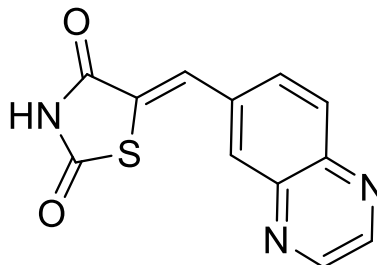
AS605240 (PI 3-Kinase Inhibitor 1)

Catalog number: B-0301

Molecular Formula: $C_{12}H_7N_3O_2S$

MW: 257.3

CAS: 648450-29-7



Alternate Name: 5-quinoxilin-6-methylene-1,3-thiazolidine-2,4-dione

Solubility: water: insoluble, DMSO: 2.5 mg/mL

Storage and Handling: Store dry between 2-8°C. Stock solutions should be stored frozen (-20°C or below).

Background: AS605240 is an ATP-competitive inhibitor of PI3-K γ (IC₅₀ = 8 nM). It is selective for the γ -isoform (IC₅₀: PI 3-K α = 60 nM, PI 3-K β = 270 nM, PI 3-K δ = 300 nM) and shows no notable activity against a wide array of protein kinases at 1 μ M. In addition, it suppresses the progression of joint inflammation in a mouse model of rheumatoid arthritis and suppresses inflammation in a mouse model of lupus.

References: 1) M. Camps *et al.* "Blockade of PI3K γ suppresses joint inflammation and damage in mouse models of rheumatoid arthritis" *Nature Medicine*, 2005, 11, 936-943. 2) D.F. Barber, *et al.*, "PI3K γ inhibition blocks glomerulonephritis and extends lifespan in a mouse model of systemic lupus" *Nature Medicine*, 2005, 11, 933-935.

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.

