



Echelon Biosciences Inc.
675 Arapen Drive, Ste 302
Salt Lake City, UT 84108
Phone: 801-588-0455
Fax: 801-588-0497
Email: echelon@echelon-inc.com
<http://www.echelon-inc.com>

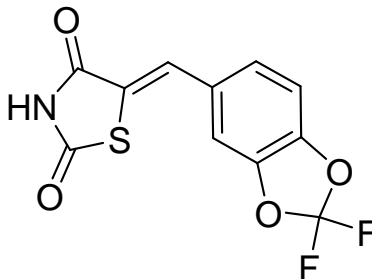
Technical Data Sheet

For research use only

Not intended or approved for
diagnostic or therapeutic use.

Product Name: AS604850 (PI 3-Kinase γ Inhibitor 2); 5-(2,2-Difluoro-benzo[1,3]dioxol-5-ylmethylene)-thiazolidine-2,4-dione

Product Number: B-0302



Formula: C₁₁H₇F₂NO₄S, FW: 287.2

Solubility: water: insoluble, DMSO: 5 mg/mL

Storage: store dry between 2-8°C. Stock solutions should be stored frozen (-20°C).

Field of Interest: 5-(2,2-Difluoro-benzo[1,3]dioxol-5-ylmethylene)-thiazolidine-2,4-dione is a selective inhibitor of PI 3-K γ with an IC₅₀ = 250 nM. It is selective for the γ -isoform (IC₅₀: PI 3-K α = 4.5 μ M, PI 3-K β , PI 3-K δ > 20 μ M) and shows no notable activity against a wide array of protein kinases at 1 μ M.

Reference: 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.

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. 5-(2,2-Difluoro-benzo[1,3]dioxol-5-ylmethylene)-thiazolidine-2,4-dione 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.