

Echelon Biosciences Inc.

LY294002

Catalog number: B-0294

Molecular Formula: C₁₉H₁₇NO₃.HCl

MW: 343.8

CAS: 154447-36-6

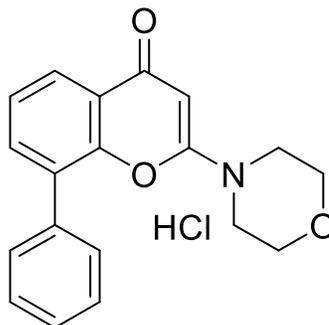
Alternate Name: 2-(4-Morpholinyl)-8-phenyl-4H-1-benzopyran-4-one

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

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

Background: LY294002 is a cell permeable, potent, and specific inhibitor of phosphatidylinositol 3-kinase (IC₅₀ = 1.4 μM; purified PI 3-kinase). It does not significantly inhibit PKC, PKA, MAP kinase, S6 kinase, EGFRK, c-src kinase, PI 4-kinase, DAG kinase or rabbit kidney ATPase at concentrations of 50 μM. Induces apoptosis. In intact neutrophils stimulated with fMet-Leu-Phe, LY294002 completely abolished PI 3-kinase activity at 50 μM without cell toxicity. The product is not sterile.

References: 1) Walker EH, Pacold ME, Perisic O, Stephens L, Hawkins PT, Wymann MP, Williams RL *Mol. Cell.* 2000, 6(4):909-19. Structural determinants of phosphoinositide 3-kinase inhibition by wortmannin, LY294002, quercetin, myricetin, and staurosporine. 2) Prestwich GD *Chemtracts - Organic Chemistry* 1994, 7, 301-305. 3)



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.

Technical Data Sheet, Rev 3, 10-12-18 – For research use only. Not intended for diagnostic or therapeutic use.



Echelon Biosciences Inc.

Ph: 866-588-0455
Fax: 801-588-0497

Echelon-inc.com