



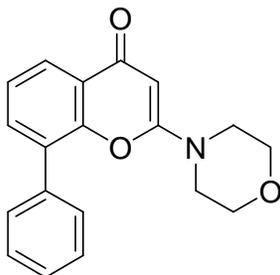
Echelon Biosciences Inc.  
675 Arapeen Drive, Suite 302  
Salt Lake City, UT 84108  
Telephone 866-588-0455  
Fax 801-588-0497  
echelon@echelon-inc.com  
www.echelon-inc.com

## Technical Data Sheet

**For research use only**  
Not intended or approved for  
diagnostic or therapeutic use.

**Product Name: 2-(4-Morpholinyl)-8-phenyl-4H-1-benzopyran-4-one (LY294002)**

**Product Number: B-0294**



**Formula:** C<sub>19</sub>H<sub>17</sub>NO<sub>3</sub>.HCl      **MW (g/mol):** 343.8

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

**Specifications and Storage Recommendations:** LY294002 should be stored dry between 2-8°C

**Field of Interest:** LY294002 is a cell permeable, potent, and specific inhibitor of phosphatidylinositol 3-kinase (IC<sub>50</sub> = 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.

**Hazardous Properties and Cautions:** The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. LY 294002 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.

### 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

**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.

Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications and methods and use, may be covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our trademarks and technologies at busdev@echelon-inc.com.