

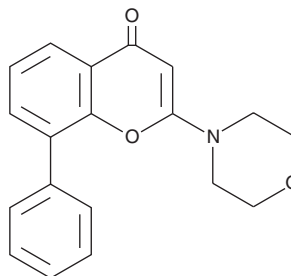
PRODUCT INFORMATION



LY294002

Item No. 70920

CAS Registry No.: 154447-36-6
Formal Name: 2-(4-morpholinyl)-8-phenyl-4H-1-benzopyran-4-one
MF: C₁₉H₁₇NO₃
FW: 307.3
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 302 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years
Melting Point: 182-184°C



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

LY294002 is supplied as a crystalline solid. LY294002 is soluble in solvents such as DMSO, ethanol, and dimethyl formamide. The solubility of LY294002 in these solvents is approximately 16 mg/ml. LY294002 is stable for at least six months in these solvents if stored at -20°C.

LY294002 is sparingly soluble in aqueous solutions (<50 µg/ml in PBS pH 7.2). For maximum solubility in aqueous buffers, LY294002 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. LY294002 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant since organic solvents may have physiological effects at low concentrations. Store aqueous solutions of LY294002 on ice and use within 12 hours of preparation.

Description

LY294002 is a phosphatidylinositol 3-kinase (PI3K) inhibitor. PI3K is an enzyme that phosphorylates the D-3 position of the inositol ring in phosphoinositides, resulting in the formation of PI(3)P, PI(3,4)P₂, and PI(3,4,5)P₃. LY294002 is a selective PI3K inhibitor with a 2.7-fold greater potency than quercetin.¹ LY294002 inhibits purified PI3K with an IC₅₀ of 1.4 µM, but does not inhibit PI4K, EGF receptor, PDGF receptor, insulin receptor, c-src, MAP kinase, S6 kinase, diacylglycerol kinase, protein kinase A, protein kinase C, and ATPase.¹ LY294002 was shown to completely abolish PI3K activity in fMet-Leu-Phe-stimulated human neutrophils, as well as, inhibit proliferation of smooth muscle cells in cultured rabbit aortic segments.¹

Reference

1. Vlahos, C.J., Matter, W.F., Hui, K.Y., *et al.* A specific inhibitor of phosphatidylinositol 3-kinase, 2-(4-morpholinyl)-8-phenyl-4H-1-benzopyran-4-one (LY294002). *J. Biol. Chem.* **269**, 5241-5248 (1994).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/29/2020

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM