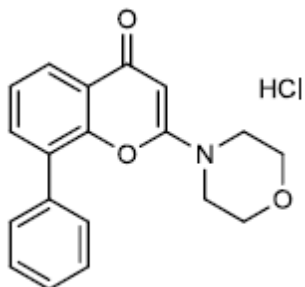


LY-294002

Cat. # AC-1005

Background: Highly selective inhibitor of phosphatidylinositol 3-kinase (IC₅₀ = 0.31, 0.73, 1.06 and 6.60 μM for PI 3-Kβ, PI 3-Kα, PI 3-Kδ and PI 3-Kγ respectively)¹. Inhibits autophagy in rat hepatocytes (IC₅₀ = 10 μM)². An extremely useful tool for determining the role of PI 3-K in cellular signaling³. Cell permeable. More stable in solution than Wortmannin, and also blocks autophagosome formation.

Application: Inhibits autophagy

**Product Information**

Purity:	>98% by TLC
MW:	307.4
Formula:	C ₁₉ H ₁₇ NO ₃
CAS No.	154447-36-6
Physical State:	Lyophilized white powder
Amount:	5mg, 25mg
Solubility:	20mg/mL in warm DMSO; 8 mg/mL in warm ethanol
Storage:	Store desiccated as supplied at -20°C for up to 3 years. Store solutions at -20°C for up to 3 months.

References

1. CJ Vlahos *et al.* *J. Biol. Chem.* 1994 269:5241
2. EF Blomaart *et al.* *Eur. J. Biochem.* 1997 243:240
3. S Chasserot-Golaz *et al.* *J. Neurochem.* 1998 70:2347

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2013 LifeSensors, Inc. All Rights Reserved