

Wortmannin

Catalog #: 27031

Lot #: 220711

Size: 10 mg

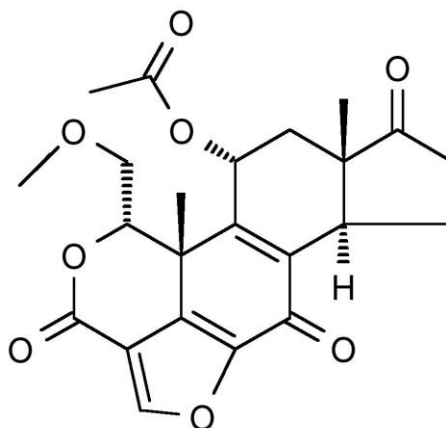
Structure:

CAS Registry #: 19545-26-7

Purity: ≥99%

Chemical Formula: C₂₃H₂₄O₈

Molecular Weight: 428.43 Da



Description: Potent and selective inhibitor of phosphatidyl-inositol 3-kinase. Active in purified preparations and cytosolic fractions (IC₅₀ = 5nM) and is highly cell permeable. Covalently binds to PI 3-kinase and is selective, inhibiting other kinases such as PI 4-kinase and myosin light chain kinase at concentrations 100-fold higher than that required for inhibition of PI 3-kinase.

Synonym(s): KY 12420

Appearance: White to off-white solid.

Solubility: Soluble in DMSO or ethanol; very poor solubility in water; maximum solubility in plain water estimated to be about 10-50µM; buffers, serum, or other additives may increase or decrease the aqueous solubility.

Storage/Stability: Store at or below -20°C. Solid form is stable at least 12 months from date of receipt, when stored as directed. Do not store aqueous solutions for more than one day. Protect from Light.

References:

1. Wymann, M.P. et al.; (1996). Mol. Cell. Biol. 16, 1722.
2. M. Ui, et al. (1995). TIBS 20, 303.
3. Arcaro, A. and Wyman, M.P. (1993). Biochem. J. 296, 297.