

Glowing products for science

Staurosporine

Staurosporine is a potent and cell-permeable inhibitor of a wide variety of protein kinases by competing for the ATP binding site.



Product attributes

Call us: 800-304-5357

CAS number	62996-74-1
Molecular weight	466.20
Storage Conditions	Store at -10 to -35 °C, Protect from light, Desiccate
Toxin	Alkaloid Toxin

Product Description

Staurosporine inhibits a wide variety of protein kinases by competing for the ATP binding site. It is a potent and cell-permeable competitive inhibitor for protein kinase C (IC50 = 0.7 nM), protein kinase A (IC50 = 7 nM), and protein kinase G (IC50 = 8.5 nM) (1,2). Staurosporine induces apoptosis in human neuroblastoma cells and chick embryonic neurons (3,4).

- Cell-permeable inhibitor of protein kinases
- Competitive inhibition at the ATP binding site
- Soluble in DMSO or ethanol
- [62996-74-1]

References

- 1. Biochem Biophys Res Commun. 158, 105 (1989), DOI: 10.1016/S0006-291X(89)80183-4
 2. Biochem Biophys Res Commun. 135, 397 (1986), DOI: 10.1016/0006-291X(86)90008-2
 3. Neuropharmacology 36, 811 (1997), DOI: 10.1046/j.1471-4159.1996.66041418.x
 4. J. Neurochem. 66, 1418 (1996), DOI: 10.1046/j.1471-4159.1996.66041418.x

This datasheet was generated on November 20, 2025 at 09:19:27 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/staurosporine/