

Staurosporine

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



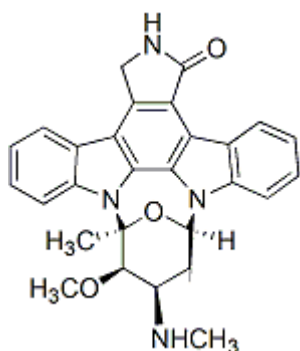
Product attributes

CAS number	62996-74-1
Molecular weight of antigen	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](https://doi.org/10.1016/S0006-291X(89)80183-4)
2. Biochem Biophys Res Commun. 135, 397 (1986), [DOI: 10.1016/0006-291X\(86\)90008-2](https://doi.org/10.1016/0006-291X(86)90008-2)
3. Neuropharmacology 36, 811 (1997), [DOI: s0028-3908\(97\)00030-0](https://doi.org/10.1016/S0028-3908(97)00030-0)
4. J. Neurochem. 66, 1418 (1996), [DOI: 10.1046/j.1471-4159.1996.66041418.x](https://doi.org/10.1046/j.1471-4159.1996.66041418.x)

This datasheet was generated on June 16, 2026 at 04:09:04 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/staurosporine/>