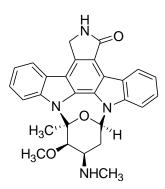


46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME:	Staurosporine
CATALOG #	00025
MOLECULAR INFORMATION:	$C_{28}H_{26}N_4O_3$ MWt: 466.53



Light yellow solid Soluble in DMSO or ethanol
Store at desiccated at -20 °C and protect from light
Staurosporine is a potent and cell-permeable inhibitor of a wide variety of protein kinases, ^{1,2} including protein kinase C ($IC_{50} = 0.7 \text{ nM}$), protein kinase A ($IC_{50} = 7 \text{ nM}$), and protein kinase G ($IC_{50} = 8.5 \text{ nM}$). Staurosporine induces apoptosis in human neuroblastoma cells and chick embryonic neurons. ^{3,4}
Ref: 1) Biochem. Biophys. Res. Commun. 158 , 105(1989); 2) Biochem. Biophys. Res. Commun. 135 , 397(1986); 3) Neuropharmacology 36 , 811(1997); 4) J. Neurochem. 66 , 1418(1996).
Not established.
Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.