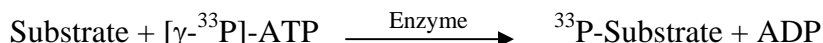


SSTK/TSSK6

(Testis-specific serine/threonine-protein kinase 6, Cancer/testis antigen 72, Serine/threonine-protein kinase SSTK, Small serine/threonine kinase, FKSG82)

- **CAT#:** SSTK/TSSK6
- **Enzyme:** Human SSTK/TSSK6
- **Substrate:** MBP (Myelin Basic Protein), 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	160
Ro 31-8220	7,400
GW 5074	65,000
H-89	>100,000
Rottlerin	>100,000

- **Scientific Information:** Q9BXA6 (<http://www.expasy.org/uniprot/Q9BXA6>)

Reference compound IC50 for SSTK/TSSK6

