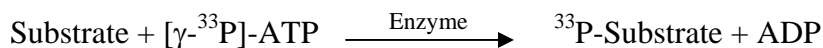


## STK22D/TSSK1

(Testis-specific serine/threonine-protein kinase 1, TSSK-1, Testis-specific kinase 1  
TSK-1, Serine/threonine-protein kinase 22A,  
TSSK1B, SPOGA1, SPOGA4, STK22A, TSSK1)

- **CAT#:** STK22D/TSSK1
- **Enzyme:** Human STK22D/TSSK1
- **Substrate:** Peptide substrate, [KKKVSRSGLYRSPSPENLNRPR], 20  $\mu$ M  
ATP 10  $\mu$ M
- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	< 0.1
Ro 31-8220	1,200
GW 5074	7,500
H-89	4,900

- **Scientific Information:** Q9BXA7 (<http://www.expasy.org/uniprot/Q9BXA7>)

### Reference Compound IC50 for STK22D

