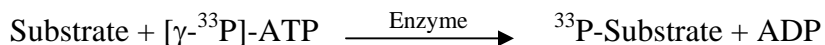


TESK1

(Dual specificity testis-specific protein kinase 1, Testicular protein kinase 1)

- **CAT#:** TESK1
- **Enzyme:** Human TESK1
- **Substrate:** Cofilin 2, 3 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	260
Ro 31-8220	22,000
GW 5074	7,500
H-89	>100,000
Rottlerin	>100,000

- **Scientific Information:** Q15569 (<http://www.expasy.org/uniprot/Q15569>)

Reference compound IC50 for TESK1

