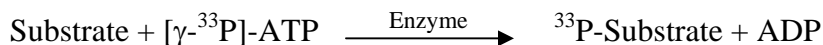


STK32C/YANK3

(Serine/threonine-protein kinase 32C, ST32C)

- **CAT#:** STK32C/YANK3
- **Enzyme:** Human STK32C/YANK3
- **Substrate:** MBP (Myelin Basic Protein), 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	250
Ro 31-8220	7,000
GW 5074	>100,000
H-89	>100,000
Rottlerin	>100,000

- **Scientific Information:** Q86UX6 (<http://www.expasy.org/uniprot/Q86UX6>)

Reference compound IC50 for STK32C/YANK3

