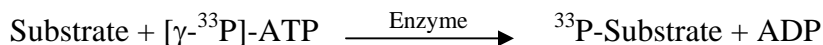


## CK2a2

(Casein kinase II subunit alpha, CSK22, CSNK2A2, CK2A2)

- **CAT#:** CK2a2
- **Enzyme:** Human CK2a2
- **Substrate:** Peptide substrate, [RRRDDDSDDD], 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	652
Ro 31-8220	72,000
GW 5074	142
Rottlerin	25,000

- **Scientific Information:** P19784 (<http://www.expasy.org/uniprot/P19784>)

### Reference Compound IC50 for CK2 $\alpha$ 2

