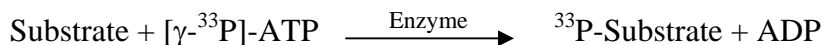


CK2a

(Casein kinase II subunit alpha, CSK21, CSNK2A1, CK2A1)

- **CAT#:** CK2a
- **Enzyme:** Human CK2a
- **Substrate:** Peptide substrate, [RRRDDDSDDD], 20 μ M
ATP 10 μ M
- **Reaction:**



- **Reference Compound IC50s:**

| Compounds | IC50 (nM) |
|---------------|-----------|
| Staurosporine | 27,600 |
| GW 5074 | 400 |
| Rottlerin | 160,000 |

- **Scientific Information:** P68400 (<http://www.expasy.org/uniprot/P68400>)

Reference compound IC50 for CK2a

