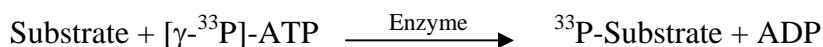


## CK1a1

(Casein kinase I isoform alpha, CKI-alpha, CK1, KC1A, CSNK1A1)

- **CAT#:** CK1a1
- **Enzyme:** Human CK1a1
- **Substrate:** Peptide substrate, [KRRRAL[pS]VASLPGL], 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	7,800
GW 5074	30,500

- **Scientific Information:** P48729 (<http://www.expasy.org/uniprot/P48729>)

### Reference Compound IC50 for CK1a1

