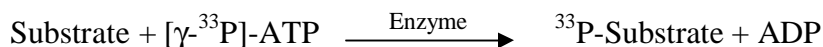


PKA

(cAMP-dependent protein kinase, PRKACA)

- **CAT#:** PKA
- **Enzyme:** Human PKA
- **Substrate:** Peptide substrate, [GRTGRRNSI-NH₂], 5 μM
ATP 1 μM

- **Reaction:**



- **Reference Compound IC₅₀s:**

Compounds	IC ₅₀ (nM)
Staurosporine	0.16
Ro 31-8220	101
GW 5074	20,000
H-89	5.5
Rottlerin	31,000

- **Scientific Information:** P17612 (<http://www.expasy.org/uniprot/P17612>)

Reference Compound IC₅₀ for PKA

