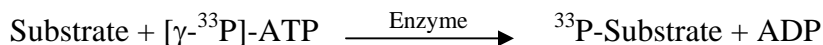


PKCb2

(Protein kinase C beta type, PKC-beta, PKC-B, PRKCB1, KPCB)

- **CAT#:** PKCb2
- **Enzyme:** Human PKCb2
- **Substrate:** Histone H1, Lipid activator, 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.8
Ro 31-8220	0.8
GW5074	17,000
H-89	2,200

- **Scientific Information:** P05771-2 (<http://www.expasy.org/uniprot/P05771-2>)

Reference compound IC50 for PKCb2

