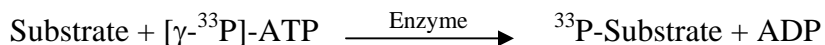


PKCd

(Protein kinase C delta type, nPKC-delta, KPCD, PRKCD)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	< 1.0
Ro 31-8220	0.65
GW5074	>20,000
H-89	280

- **Scientific Information:** Q05655 (<http://www.expasy.org/uniprot/Q05655>)

Reference compound IC50 for PKCd

