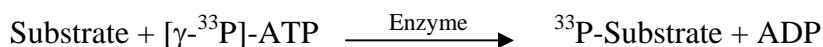


PKCtheta

(Protein kinase C theta type, nPKC-theta, PRKCQ, PRKCT, KPCT)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3.1
Ro 31-8220	3.8

- **Scientific Information:** Q04759 (<http://www.expasy.org/uniprot/Q04759>)

Reference compound IC50 for PKCtheta

