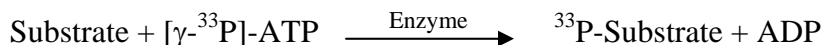


PKC ι

(Protein kinase C iota type, nPKC-iota, Atypical protein kinase C-lambda/iota, aPKC-lambda/iota, PRKC-lambda/iota, PRKCI, PKCI)

- **CAT#:** PKCiota
- **Enzyme:** Human PKCiota
- **Substrate:** Peptide substrate, [ERM RPRKRQGSVRRRV], 20 μM
ATP 1 μM

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	4.0
Ro 31-8220	76.8
H-89	11,800

- **Scientific Information:** P41743 (<http://www.expasy.org/uniprot/P41743>)

