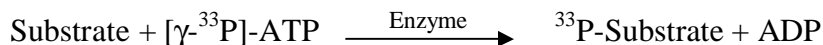


PKC μ

(Serine/threonine-protein kinase D1, nPKC-D1, Protein kinase D
Protein kinase C mu type, nPKC-mu, PRKD1, PKD, PKD1, PRKCM, KPCD1)

- **CAT#:** PKC μ
- **Enzyme:** Human PKC μ
- **Substrate:** Peptide substrate, [KKLNRTLSVA], 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.3
Ro 31-8220	62
GW 5074	6,600
H-89	562

- **Scientific Information:** Q15139 (<http://www.expasy.org/uniprot/Q15139>)

Reference Compound IC50 for PKC μ

