

## PKCmu/PRKD1

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

- **CAT#:** PKCmu
- **Enzyme:** Human PKCmu
- **Substrate:** Peptide substrate, [KKLNRTLSVA], 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.5
Ro 31-8220	210
GW 5074	5,700
H-89	910

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

### Reference compound IC50 for PKCmu/PRKD1

