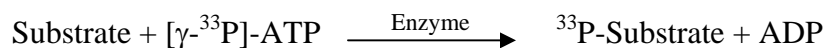


## CAMK2d

(Calcium/calmodulin-dependent protein kinase type II delta chain, CaM-kinase II delta chain  
CaM kinase II subunit delta, CaMK-II subunit delta, CAMK2D, CAMKD)

- **CAT#:** CAMK2d
- **Enzyme:** Human CAMK2d
- **Substrate:** Peptide substrate, [KKLNRTLSEFAEPG], 20  $\mu$ M  
Ca<sup>2+</sup>-Calmodulin, 1  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	< 0.5
Ro 31-8220	650
GW 5074	3,100
H-89	15,000

- **Scientific Information:** Q13557 (<http://www.expasy.org/uniprot/Q13557>)

### Reference compound IC50 for CAMK2d

