

CAMK1d

(Calcium/calmodulin-dependent protein kinase type 1D, CaM kinase ID, CaM kinase I delta, CaMKI-delta, CaM-KI delta, CaMKI delta, Camk1D, CamKI-like protein kinase, CKLiK, CAMK1D, KCC1D)

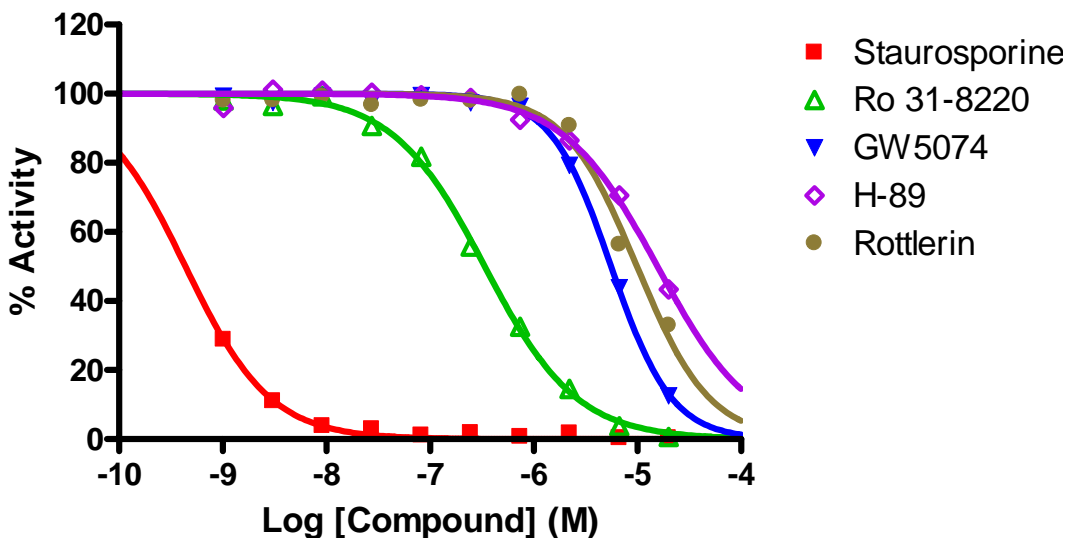
- **CAT#:** CAMK1d
- **Enzyme:** Human CAMK1d
- **Substrate:** Peptide substrate, [KKALRRQETVDAL], 20 μ M
Ca²⁺-Calmodulin, 1 μ M
ATP 10 μ M
- **Reaction:**

$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} \text{}^{33}\text{P}\text{-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.4
Ro 31-8220	340
GW 5074	5,600
H-89	16,000
Rottlerin	10,000

- **Scientific Information:** Q8IU85 (<http://www.expasy.org/uniprot/Q8IU85>)

Reference compound IC50 for CAMK1d



Service performed using kinases under license from  invitrogen™ part of *Life* technologies™