

## CAMK4

(Calcium/calmodulin-dependent protein kinase type IV, CAM kinase-GR, CaMK IV, CAMK4, KCC4)

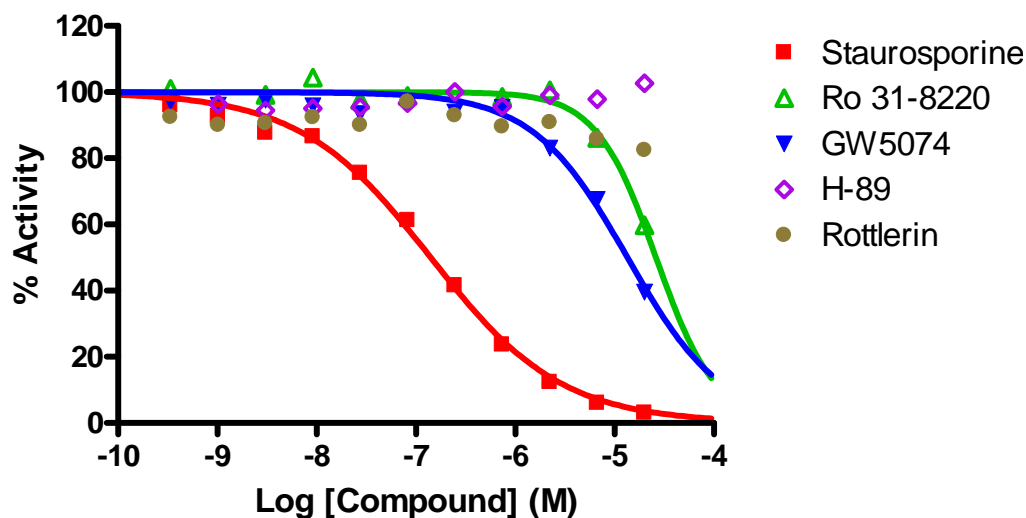
- **CAT#:** CAMK4
- **Enzyme:** Human CAMK4
- **Substrate:** Peptide substrate, [KKLNRTLSEFAEPG], 20  $\mu$ M  
Ca<sup>2+</sup>-Calmodulin, 1  $\mu$ M  
ATP 10  $\mu$ M
- **Reaction:**  

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

Compounds	IC50 (nM)
Staurosporine	141
Ro 31-8220	26,100
GW 5074	13,500

- **Scientific Information:** Q9UQM7 (<http://www.expasy.org/uniprot/Q9UQM7>)

### Reference Compound IC50 for CAMK4



Service performed  
using kinases  
under license from  **invitrogen**<sup>™</sup>  
part of *life* technologies<sup>™</sup>