

## DCAMKL2

(Doublecortin-like and CAM kinase-like 2, Doublecortin-like kinase 2, DCLK2, DCDC3B, DCK2)

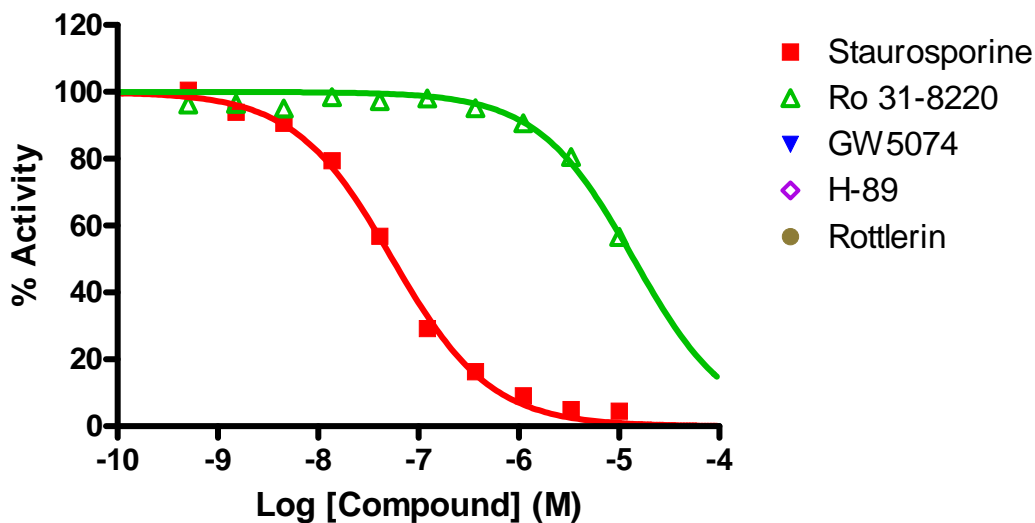
- **CAT#:** DCAMKL2
- **Enzyme:** Human DCAMKL2
- **Substrate:** Peptide substrate, [KKALRRQETVDAL], 20  $\mu$ M  
Ca<sup>2+</sup>-Calmodulin, 1  $\mu$ M  
ATP 10  $\mu$ M
- **Reaction:**  

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

Compounds	IC50 (nM)
Staurosporine	55.1
Ro 31-8220	14,000

- **Scientific Information:** Q8N568 (<http://www.expasy.org/uniprot/Q8N568>)

### Reference Compound IC50 for DCAMKL2



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