

## RIPK5

(Receptor-interacting serine/threonine-protein kinase 5, Dusty protein kinase, Dusty PK, RIP-homologous kinase, Sugen kinase 496, KIAA0472, RIP5, SGK496, HDCMD38P)

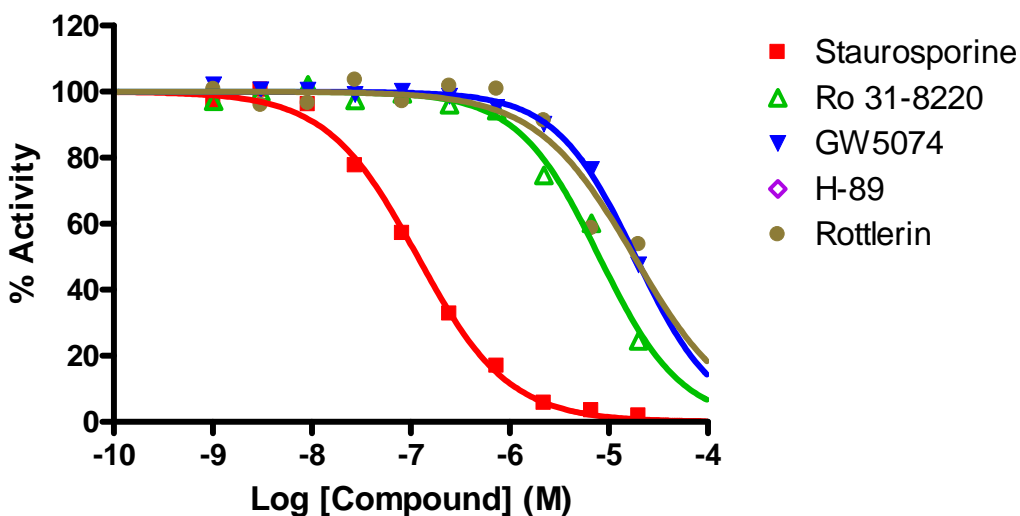
- **CAT#:** RIPK5
- **Enzyme:** Human RIPK5
- **Substrate:** MBP (Myelin Basic Protein), 10  $\mu$ M  
ATP 10  $\mu$ M
- **Reaction:**  

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

Compounds	IC50 (nM)
Staurosporine	120
Ro 31-8220	8,000
GW 5074	19,000
Rottlerin	20,000

- **Scientific Information:** Q6XUX3 (<http://www.expasy.org/uniprot/Q6XUX3>)

### Reference Compound IC50 for RIPK5



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