

## MNK2

(MAP kinase-interacting serine/threonine-protein kinase 2,  
MAP kinase signal-integrating kinase 2, Mnk2, MKNK2)

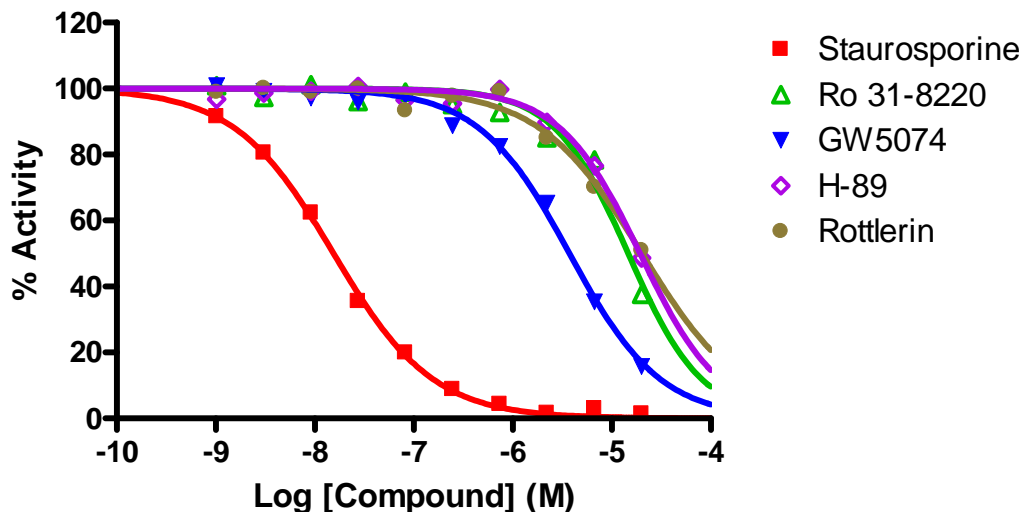
- **CAT#:** MNK2
- **Enzyme:** Human MNK2
- **Substrate:** MBP (Myelin Basic Protein), 10  $\mu$ M  
ATP 10  $\mu$ M
- **Reaction:**  

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

Compounds	IC50 (nM)
Staurosporine	15.5
Ro 31-8220	14,600
GW 5074	3,700
H-89	19,300
Rottlerin	20,000

- **Scientific Information:** Q9HBH9 (<http://www.expasy.org/uniprot/Q9HBH9>)

### Reference compound IC50 for MNK2/MKNK2



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