

## MLCK

(Myosin light chain kinase-smooth muscle, Telokin, Kinase-related protein, KRP, MYLK)

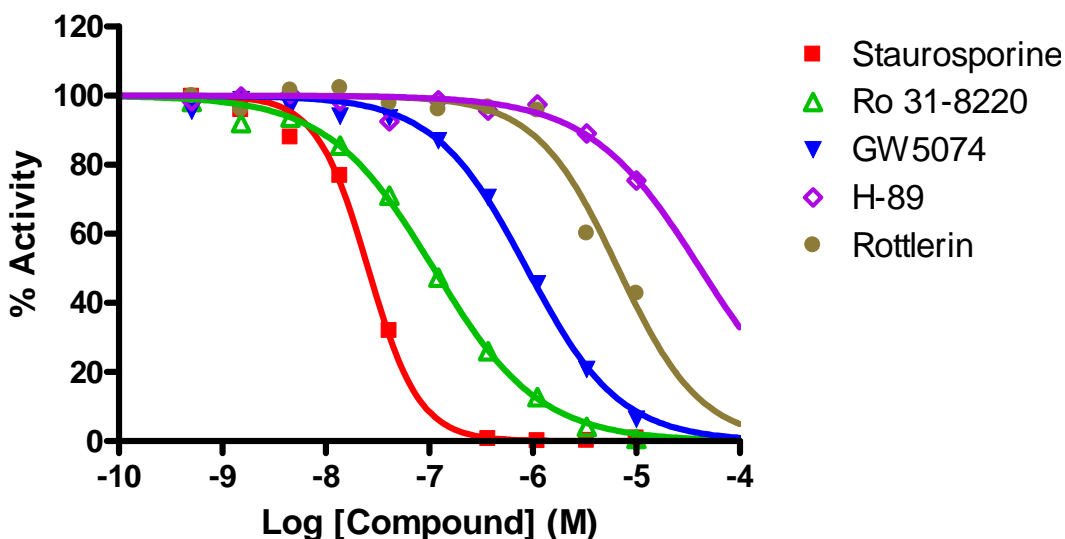
- **CAT#:** MLCK
- **Enzyme:** Human MLCK
- **Substrate:** Peptide substrate, [KKLNRTLSFAEPG], 20  $\mu$ M  
Ca<sup>2+</sup>-Calmodulin, 1  $\mu$ M  
ATP 10  $\mu$ M
- **Reaction:**  

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

| Compounds     | IC50 (nM) |
|---------------|-----------|
| Staurosporine | 25.6      |
| Ro 31-8220    | 110       |
| GW 5074       | 876       |
| H-89          | 41,600    |
| Rottlerin     | 7,000     |

- **Scientific Information:** Q15746 (<http://www.expasy.org/uniprot/Q15746>)

### Reference compound IC50 for MLCK/MYLK



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