

## ERK5/MAPK7

(Mitogen-activated protein kinase 7, Big MAP kinase 1, Extracellular signal-regulated kinase 5, BMK1, PRKM7, MK07)

- **CAT#:** ERK5/MAPK7
- **Enzyme:** Human ERK5/MAPK7
- **Substrate:** MBP (Myelin Basic Protein), 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	122
Ro 31-8220	66
GW5074	310
H-89	>50,000
Rottlerin	>50,000

- **Scientific Information:** Q13164 (<http://www.expasy.org/uniprot/Q13164>)

### Reference compound IC50 for ERK5/MAPK7

