

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 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	10,000
Ro 31-8220	49,000
GW5074	9,100
H-89	>50,000
Rottlerin	>50,000

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

Reference compound IC50 for ERK5/MAPK7

