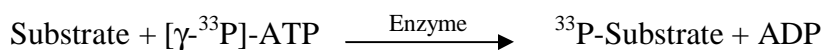


## ERK1

(Mitogen-activated protein kinase 3, Extracellular signal-regulated kinase 1, Insulin-stimulated MAP2 kinase, MAP kinase 1, MAPK 1, p44-ERK1, ERT2, p44-MAPK, Microtubule-associated protein 2 kinase, MK03)

- **CAT#:** ERK1
- **Enzyme:** Human ERK1
- **Substrate:** MBP (Myelin Basic Protein), 5  $\mu$ M  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	5,100
Ro 31-8220	19,000
Rottlerin	9,200

- **Scientific Information:** P27361 (<http://www.expasy.org/uniprot/P27361>)

