

ERK2

(Mitogen-activated protein kinase 2, Extracellular signal-regulated kinase 2
MAP kinase 2, p42-MAPK, ERT1, MK01)

- **CAT#:** ERK2
- **Enzyme:** Human ERK2
- **Substrate:** MBP (Myelin Basic Protein), 5 μ M
ATP 1 μ M
- **Reaction:**
$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} {}^{33}\text{P}\text{-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	522
Ro 31-8220	18,500
GW 5074	14,700
H-89	36,400
Rottlerin	6,800

- **Scientific Information:** P28482 (<http://www.expasy.org/uniprot/P28482>)

Reference compound IC50 for ERK2

