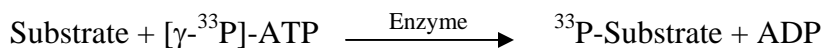


P38g/MAPK12

(Mitogen-activated protein kinase 12, Mitogen-activated protein kinase p38 gamma, Extracellular signal-regulated kinase 6, ERK-6, ERK5, Stress-activated protein kinase 3, MAP kinase p38 gamma, MK12)

- **CAT#:** P38g/MAPK12
- **Enzyme:** Human P38g/MAPK12
- **Substrate:** MBP (Myelin Basic Protein), 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	4,400
GW 5074	>20,000
Rottlerin	>20,000

- **Scientific Information:** P53778 (<http://www.expasy.org/uniprot/P53778>)

Reference compound IC50 for P38g/MAPK12

