

---

## p38γ(MAPK12)

Product code 04-155

Full-length human p38γ [1-367(end) amino acids of accession number NP\_002960.2] was expressed as N-terminal GST-fusion protein (69 kDa) using E. coli expression system. GST-p38γ was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K6. Activated GST-p38γ was purified using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Modified Erktide

ATP (μM) Km app / Bin : 13 / 10

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 88

IC50 at 1 mM ATP (nM) : 2800

---