

---

---

## p38 $\gamma$ (MAPK12)

Product code: 04-155

Full-length human p38 $\gamma$  [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-p38 $\gamma$  was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K6. Activated GST-p38 $\gamma$  was purified using glutathione sepharose chromatography.

Assay platform	: MSA
Substrate	: Modified Erktide
ATP ( $\mu$ M) Km app / Bin	: 13 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 88
IC50 at 1 mM ATP (nM)	: 2800

---

---