

---

---

## p38 $\delta$ (MAPK13)

Product code: 04-154

Full-length human p38 $\delta$  [1-365(end) amino acids of accession number NP\_002745.1] was expressed as N-terminal GST-fusion protein (69 kDa) using E. coli expression system. GST-p38 $\delta$  was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K6. Activated GST-p38 $\delta$  was purified using glutathione sepharose chromatography.

Assay platform	: MSA
Substrate	: Modified Erktide
ATP ( $\mu$ M) Km app / Bin	: 5.8 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 220
IC50 at 1 mM ATP (nM)	: >10000

---

---