
p38δ(MAPK13)

Product code 04-154

Full-length human p38δ [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-p38δ was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K6. Activated GST-p38δ was purified using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Modified Erktide

ATP (μM) Km app / Bin : 5.8 / 5

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 220

IC50 at 1 mM ATP (nM) : >10000
