
p38 β (MAPK11)

Product code 04-153

Full-length human p38 β [1-364(end) amino acids of accession number NP_002742.3] was expressed as N-terminal GST-fusion protein (69 kDa) using E. coli expression system. GST-p38 β was purified by using glutathione chromatography.

Assay platform : Mobility Shift Assay

Substrate : Modified Erktide

ATP (μ M) Km app / Bin : 63 / 75

Metal : Mg

Reference compound : SB202190

IC50 at ATP Bin (nM) : 16

IC50 at 1 mM ATP (nM) : 110
