
JNK3(MAPK10)

Product code 04-150

Full-length human JNK3 [1-426(end) amino acids of accession number NP_620446.1] was expressed as N-terminal GST-fusion protein (75 kDa) using E.coli expression system. GST-JNK3 was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K4 and MAP2K7. Activated GST-JNK3 was purified using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Modified Erktide

ATP (μ M) Kmapp / Bin : 6 / 25

Metal : Mg

Reference compound : K252a

IC50 at ATP Bin (nM) : 26

IC50 at 1 mM ATP (nM) : 730
