
JNK2(MAPK9)

Product code 04-164

Human JNK2, catalytic domain [1-364 amino acids of accession number NP_002743.3] was expressed as N-terminal GST-fusion protein (69 kDa) using E. coli expression system. GST-JNK2 was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K4 and MAP2K7. Activated GST-JNK2 was purified using Ni-NTA chromatography.

Assay platform : Mobility Shift Assay

Substrate : Modified Erktide

ATP (μ M) K_{mapp} / B_{in} : 21 / 50

Metal : Mg

Reference compound : K252a

IC₅₀ at ATP B_{in} (nM) : 110

IC₅₀ at 1 mM ATP (nM) : 1600
