
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	: MSA
Substrate	: Modified Erktide
ATP (μ M) Km app / Bin	: 6 / 25
Metal	: Mg
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 26
IC50 at 1 mM ATP (nM)	: 730
