

---

## 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	: MSA
Substrate	: Modified Erktide
ATP ( $\mu$ M) Km app / Bin	: 21 / 50
Metal	: Mg
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 110
IC50 at 1 mM ATP (nM)	: 1600

