

---

# JNK1(MAPK8)

Product code: 04-163

Human JNK1, catalytic domain [2-364 amino acids of accession number NP\_620634.1] was expressed as N-terminal GST-fusion protein (69 kDa) using E. coli expression system. GST-JNK1 was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K4 and MAP2K7. Activated GST-JNK1 was purified using glutathione sepharose chromatography.

Assay platform : MSA  
Substrate : Modified Erktide

ATP ( $\mu$ M) Km app / Bin : 29 / 100  
Metal : Mg  
Reference compound : K252a  
IC50 at ATP Bin (nM) : 99  
IC50 at 1 mM ATP (nM) : 770

---