
p38 α (MAPK14)

Product code: 04-152

Truncated human p38 α [9-352 amino acids of accession number NP_620581.1] was expressed as N-terminal GST-fusion protein (66 kDa) using E. coli expression system. GST-p38 α was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K6. Activated GST-p38 α was purified using glutathione sepharose chromatography.

Assay platform	: MSA
Substrate	: Modified Erktide
ATP (μ M) Km app / Bin	: 150 / 150
Metal	: Mg
Reference compound	: SB202190
IC50 at ATP Bin (nM)	: 6.3
IC50 at 1 mM ATP (nM)	: 22
