
Erk2(MAPK1)

Product code: 04-143

Full-length human Erk2 [1-360(end) amino acids of accession number NP_002736.3] was expressed as N-terminal GST-fusion protein (69 kDa) using E.coli expression system. GST-Erk2 was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K1. Activated GST-Erk2 was purified using glutathione sepharose chromatography.

Assay platform : MSA
Substrate : Modified Erktide

ATP (μ M) Km app / Bin : 33 / 50
Metal : Mg
Reference compound : K252a
IC50 at ATP Bin (nM) : 21
IC50 at 1 mM ATP (nM) : 180
