
Erk1(MAPK3)

Product code 04-142

Full-length human Erk1 [1-379(end) amino acids of accession number NP_002737.1] was expressed as N-terminal GST-fusion protein (70 kDa) using E.coli expression system. GST-Erk1 was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K1. Activated GST-Erk1 was purified using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Modified Erktide

ATP (μ M) Kmapp / Bin : 34 / 50

Metal : Mg

Reference compound : K252a

IC50 at ATP Bin (nM) : 37

IC50 at 1 mM ATP (nM) : 400
