

---

## Erk1(MAPK3)

Product code: 04-142

Full-length human Erk1 [1-379(end) amino acids of accession number NP\_002737.2] 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 : MSA  
Substrate : Modified Erktide

ATP ( $\mu$ M) Km app / Bin : 34 / 50

Metal : Mg

Reference compound : K252a

IC50 at ATP Bin (nM) : 37

IC50 at 1 mM ATP (nM) : 400

---