

---

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

---