

---

## Erk5(MAPK7)

Product code: 04-146

Human Erk5, catalytic domain [1-398 amino acids of accession number NP\_002740.2] was expressed as N-terminal GST-fusion protein (72 kDa) using E. coli expression system. GST-Erk5 was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K5. Activated GST-Erk5 was purified by using Ni-NTA affinity chromatography.

Assay platform	: MSA
Substrate	: EGFR-derived peptide
ATP ( $\mu$ M) Km app / Bin	: 450 / 1000
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 280
IC50 at 1 mM ATP (nM)	: 280

---