

---

# MAP2K2

Product code: 07-142

Full-length human MAP2K2 [1-400(end) amino acids of accession number NP\_109587.1] was co-expressed as N-terminal GST-fusion protein (71 kDa) with human His-tagged RAF1 [306-648(end) amino acids and Y340D and Y341D of accession number NP\_002871.1] using baculovirus expression system.

Assay platform	:	MSA
Substrate	:	Erk2
		Cascade Assay *
ATP ( $\mu$ M)	:	1000
Metal	:	Mg
Reference compound	:	Staurosporine
IC50 at 1 mM ATP (nM)	:	54

\* Erk2/Modified Erktide

---