

---

# 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. GST-MAP2K2 was purified by using glutathione sepharose chromatography and Ni-NTA affinity chromatography.

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

---