

---

---

# MAP2K5

Product code: 07-145

Full-length human MAP2K5 [1-448(end) amino acids of accession number NP\_660143.1] was co-expressed as N-terminal GST-fusion protein (77 kDa) with human His-tagged MAP3K3[1-626(end) amino acids of accession number NP\_002392.2] , CDC37 and HSP90 using baculovirus expression system.

Assay platform	: MSA
Substrate	: Erk5
	Cascade Assay *
ATP ( $\mu$ M)	: 1000
Metal	: Mg
Reference compound	: Staurosporine
IC50 at 1 mM ATP (nM)	: 62

\* Erk5/EGFR-derived peptide

---

---