

---

# MAP4K2

Product code: 07-111

Full-length human MAP4K2 [1-820(end) amino acid of accession number NP\_004570.2] was expressed as N-terminal GST-fusion protein (119 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: S6K2 peptide
ATP ( $\mu$ M) Km app / Bin	: 93 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.3
IC50 at 1 mM ATP (nM)	: 0.87

