

---

---

## HGK(MAP4K4)

Product code: 07-137

Human HGK, catalytic domain [1-328 amino acids of accession number NP\_004825.2] was expressed as N-terminal GST-fusion protein (64 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Moesin-derived peptide
ATP ( $\mu$ M) Km app / Bin	: 9.4 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.0
IC50 at 1 mM ATP (nM)	: 23

---

---