

---

## ***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. GST-HGK was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay  
Substrate : Moesin-derived peptide

ATP ( $\mu$ M)  $K_{mapp}$  /  $B_{in}$  : 9.4 / 10

Metal : Mg

Reference compound : Staurosporine

IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 1.0

IC<sub>50</sub> at 1 mM ATP (nM) : 23

---