

---

# HIPK1

Product code 04-135

Human HIPK1, catalytic domain [158-555 amino acids of accession number NP\_689909.2] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-HIPK1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay  
Substrate : DYRKtide-F

ATP ( $\mu$ M) Km app / Bin : 4.4 / 5

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 570

IC50 at 1 mM ATP (nM) : n.a.

---