

---

---

# HIPK3

Product code: 04-137

Human HIPK3, catalytic domain [161-562 amino acids of accession number NP\_005725.3] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: DYRKtide-F
ATP ( $\mu$ M) Km app / Bin	: 7.3 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 120
IC50 at 1 mM ATP (nM)	: >10000

