

---

---

# HIPK2

Product code: 04-136

Full-length human HIPK2 [1-1198(end) amino acids of accession number NP\_073577.3] was expressed as N-terminal GST-fusion protein (158 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: DYRKtide-F
ATP ( $\mu$ M) Km app / Bin	: 5.9 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 170
IC50 at 1 mM ATP (nM)	: n.a.

