
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. GST-HIPK2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : DYRKtide-F

ATP (μ M) K_m app / B_{in} : 5.9 / 5

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 170

IC₅₀ at 1 mM ATP (nM) : n.a.
