
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. GST-HIPK3 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : DYRKtide-F

ATP (μ M) K_m app / B_{in} : 7.3 / 5

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 120

IC₅₀ at 1 mM ATP (nM) : >10000
