
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 (μ M) K_m app / Bin : 7.3 / 5
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 120
IC50 at 1 mM ATP (nM) : >10000
