
HIPK4

Product code: 04-138

Full-length human HIPK4 [1-616(end) amino acids of accession number NP_653286.2] was expressed as N-terminal GST-fusion protein (96 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : DYRKtide-F

ATP (μ M) Km app / Bin : 7 / 5

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 71

IC50 at 1 mM ATP (nM) : >10000
