
HIPK1

Product code: 04-135

Human HIPK1, catalytic domain [158-555 amino acids of accession number NP_689909.2] 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 / B_{in} : 4.4 / 5
Metal : Mg
Reference compound : Staurosporine
IC₅₀ at ATP B_{in} (nM) : 570
IC₅₀ at 1 mM ATP (nM) : n.a.
