
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.

