
PHKG1

Product code: 02-152

Full-length human PHKG1 [1-387(end) amino acids of accession number NP_006204.1] was expressed as N-terminal GST-fusion protein (72 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: GS peptide
ATP (μ M) Km app / Bin	: 71 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.22
IC50 at 1 mM ATP (nM)	: n.a.

