
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.
