
PHKG2

Product code: 02-153

Full-length human PHKG2 [1-406(end) amino acids of accession number NP_000285.1] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: GS peptide
ATP (μ M) Km app / Bin	: 8.1 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.74
IC50 at 1 mM ATP (nM)	: n.a.

