
PDK1(PDPK1)

Product code: 01-132

Full-length human PDK1 [1-556(end) amino acids of accession number NP_002604.1] was expressed as N-terminal GST-fusion protein (91 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: T308tide
ATP (μ M) Km app / Bin	: 9.6 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 9.2
IC50 at 1 mM ATP (nM)	: 12

