
PAK2

Product code: 07-124

Full-length human PAK2 [1-524(end) amino acids of accession number NP_002568.2] was expressed as N-terminal GST-fusion protein (85 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : DAPK1tide

ATP (μ M) Km app / Bin : 81 / 100
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 4.5
IC50 at 1 mM ATP (nM) : 22
