
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. GST-PAK2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : DAPK1tide

ATP (μ M) K_{mapp} / B_{in} : 81 / 100
Metal : Mg
Reference compound : Staurosporine
IC₅₀ at ATP B_{in} (nM) : 4.5
IC₅₀ at 1 mM ATP (nM) : 22
