
PAK5(PAK7)

Product code 07-127

Human PAK5, catalytic domain [425-719(end) amino acids of accession number NP_065074.1] was expressed as N-terminal GST-fusion protein (60 kDa) using baculovirus expression system. GST-PAK5 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : DAPK1tide

ATP (μ M) Km app / Bin : 1.9 / 1

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 2.5

IC50 at 1 mM ATP (nM) : 290
