
PAK1

Product code 07-123

Full-length human PAK1 [1-545(end) amino acids of accession number NP_002567.3] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-PAK1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : LIMKtide

ATP (μ M) Km app / Bin : 300 / 300

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 4.0

IC50 at 1 mM ATP (nM) : 11
