
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.

Assay platform : MSA
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
