
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
