
PAK6

Product code: 07-128

Full-length human PAK6 [1-681(end) amino acids of accession number NP_064553.1] was expressed as N-terminal GST-fusion protein (102 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: SGKtide
ATP (μ M) Km app / Bin	: 3.7 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.2
IC50 at 1 mM ATP (nM)	: n.a.

