

---

---

## 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.

Assay platform	: MSA
Substrate	: DAPK1tide
ATP ( $\mu$ 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

