

---

# PAK2

Product code: 07-124

Full-length human PAK2 [1-524(end) amino acids of accession number NP\_002568.2] was expressed as N-terminal GST-fusion protein (85 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : DAPK1tide  
  
ATP ( $\mu$ M)  $K_m$  app / Bin : 81 / 100  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 4.5  
IC50 at 1 mM ATP (nM) : 22

---