

---

# 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 ( $\mu$ 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

---