

---

# TSSK1

Product code: 02-364

Full-length human TSSK1 [1-367(end) amino acids of accession number NP\_114417.1] was expressed as N-terminal GST-fusion protein using baculovirus expression system. GST-TSSK1 was purified by using glutathione sepharose chromatography. GST-TSSK1 was cleaved by PreScission protease and GST-free TSSK1 (42 kDa) was collected as flow-through fraction from glutathione sepharose chromatography.

Assay platform : MSA  
Substrate : GS peptide  
  
ATP ( $\mu$ M) Km app / Bin : 11 / 10  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 0.19  
IC50 at 1 mM ATP (nM) : 0.95

---