
TSSK1B(TSSK1)

Product code: 02-364

Full-length human TSSK1B [1-367(end) amino acids of accession number NP_114417.1] was expressed as N-terminal GST-fusion protein using baculovirus expression system. GST-TSSK1B was purified by using glutathione sepharose chromatography. GST-TSSK1B was cleaved by PreScission protease and GST-free TSSK1B (42 kDa) was collected as flow-through fraction from glutathione sepharose chromatography.

Assay platform : MSA
Substrate : GS peptide

ATP (μ 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
