
TSSK2

Product code: 02-165

Full-length human TSSK2 [1-358(end) amino acids of accession number NP_443732.3] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : GS peptide

ATP (μ M) Km app / Bin : 8.8 / 10
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 4.7
IC50 at 1 mM ATP (nM) : n.a.
