

---

# 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 ( $\mu$ 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.

