
TSSK3

Product code: 02-166

Full-length human TSSK3 [2-268(end) amino acids of accession number NP_443073.1] was expressed as N-terminal GST-fusion protein (57 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: GS peptide
ATP (μ M) Km app / Bin	: 45 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 12
IC50 at 1 mM ATP (nM)	: n.a.

