
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.
