
AurA(AURKA)/TPX2

Product code: 05-101

Full-length human AurA [1-403(end) amino acids of accession number NP_940835.1] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : Kemptide

ATP (μ M) Km app / Bin : 1.7 / 2
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 6.1
IC50 at 1 mM ATP (nM) : n.a.
