
AurA(AURKA)/TPX2

Product code 05-186

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. GST-AurA was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Kemptide

ATP (μ M) K_{mapp} / B_{in} : 1.7 / 2

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 6.1

IC₅₀ at 1 mM ATP (nM) : n.a.
