

---

---

## 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 ( $\mu$ 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.

