

---

## AurA(AURKA)

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 : 27 / 25  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 1.8  
IC50 at 1 mM ATP (nM) : 17

---