

---

## AurB(AURKB)/INCENP

Product code: 05-102

Full-length human AurB [1-344(end) amino acids and M298T of accession number NP\_004208.2] was co-expressed as N-terminal GST-fusion protein (66 kDa) with His-tagged INCENP(INBOX) [803-918(end) amino acids of accession number NP\_001035784.1] using baculovirus expression system. GST-AurB was purified by using glutathione sepharose chromatography.

Assay platform : MSA  
Substrate : Kemptide

ATP ( $\mu$ M) Km app / Bin : 16 / 25  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 7.1  
IC50 at 1 mM ATP (nM) : 62

---