

---

---

# 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.

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

