

Product Information

AurB(AURKB)/INCENP

Product Number : 05-102

Product description

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.

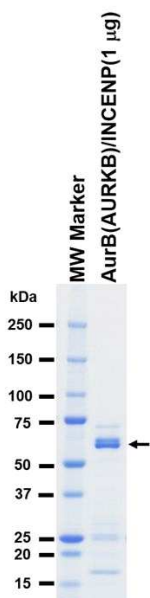
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

Store at -80C.
Avoid repeating freeze-thaws.

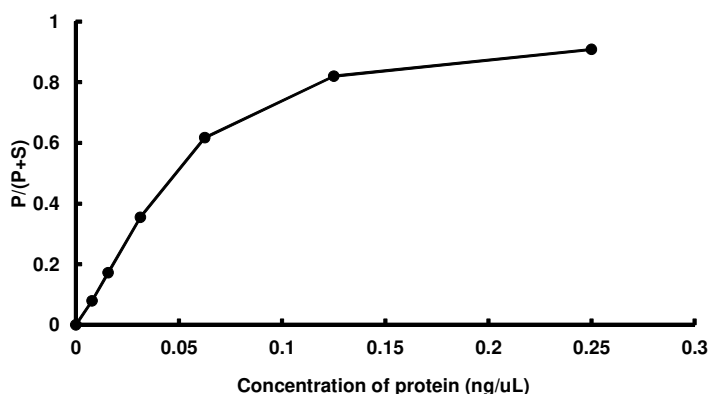
SDS-PAGE



Purity: 68 %

The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was measured by off-chip mobility shift assay(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : Kemptide

ATP : 100 μM