

## Product Information

### BTN-AurB(AURKB)/INCENP

 Product Number : **05-402-21N**

#### Product description

Full-length human AurB [1-344(end) amino acids of accession number Q96GD4] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (43 kDa) with His-tagged INCENP(INBOX) [803-918(end) amino acids of accession number AAU04398] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose and activated with ATP. Activated protein was purified by using gel filtration.

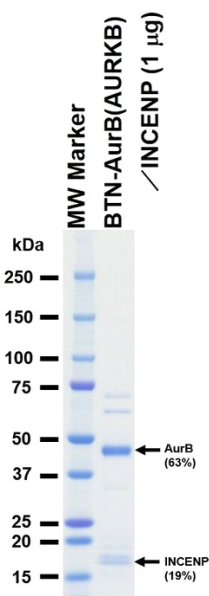
#### Storage buffer:

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.

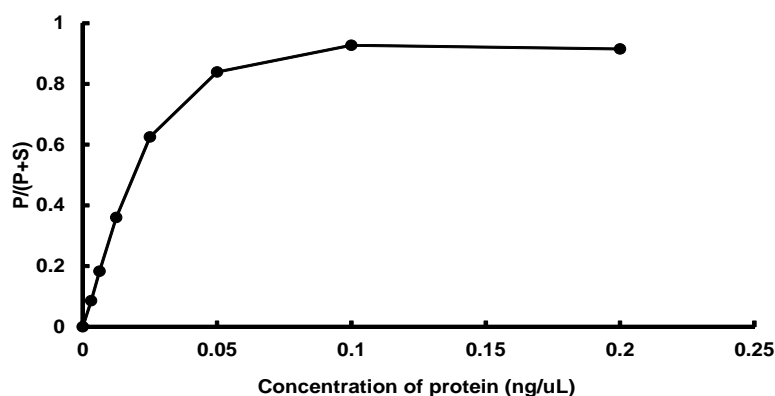
#### SDS-PAGE



Purity: 82 %

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  $\mu$ M