

## Product Information

### AurA(AURKA)

 Product Number : **05-101**

#### Product description

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. GST-AurA was purified by using glutathione sepharose chromatography.

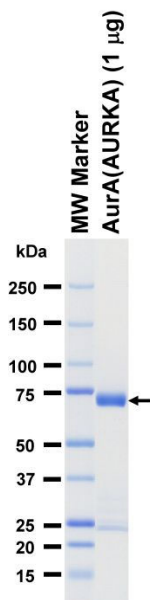
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,  
 1 mM DTT, 10% glycerol, pH 7.5

#### Storage and Handling:

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

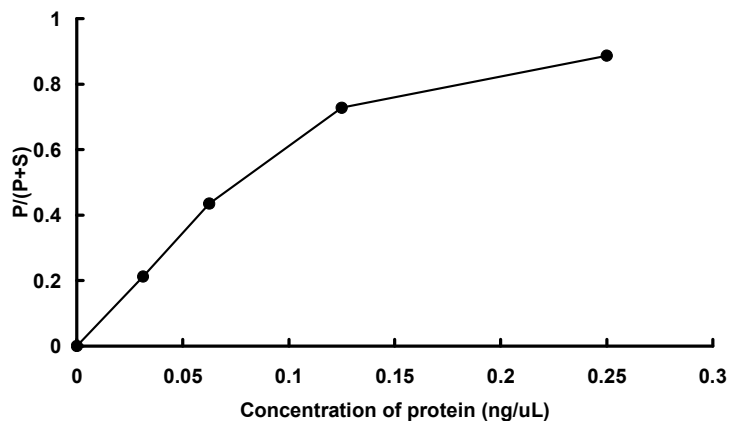
#### SDS-PAGE



Purity: 86 %

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

#### Activity data



The activity was measured by off-chip mobility shift assay. The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by LabChip™ 3000.

Substrate : Kemptide

ATP : 100 µM