

## AURORA B, Unactive

Full-length recombinant protein expressed in Sf9 cells

**Catalog # A31-14G**

Lot # X635-2

### Product Description

Recombinant full-length human AURORA B was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_004217](#). It can be activated by INCENP *in vitro*.

### Gene Aliases

AURKB, AIK2; AIM1; ARK2; AurB; IPL1; AIM-1; STK12

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

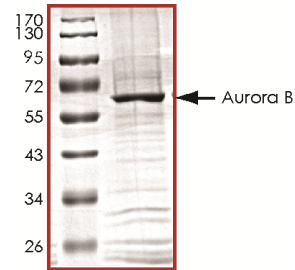
### Scientific Background

AURORA B is a member of the Aurora kinase family that associates with microtubules during chromosome movement and segregation. AURORA B localizes to the microtubules near kinetochores, specifically to the specialized microtubules called K-fibers (1). AURORA B inhibits the microtubule depolymerizing activity of mitotic centromere-associated kinesin (MCAK) by phosphorylating MCAK on Ser92 (2). This phosphorylation also regulates MCAK translocalization from kinetochores to the centromere. AURORA B has been identified as a target for the development of new anticancer agents since inhibition of AURORA B gives rise to the more pronounced antiproliferative phenotype.

### References

- Shannon, K B. et al: Chromosome dynamics: new light on Aurora B kinase function. *Curr Biol.* 2002 Jul 9;12(13):R458-60.
- Andrews, P. D. et al: Aurora B regulates MCAK at the mitotic centromere. *Dev. Cell* 6: 253-268, 2004.

### Purity



The purity of Aurora B was determined to be **>75%** by densitometry. Approx. MW **68kDa**.

## AURORA B, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog Number **A31-14G**

Specific Lot Number **X635-2**

Purity **>75%**

Concentration **0.1 µg/µl**

Stability **1yr At  $-70^{\circ}\text{C}$  from date of shipment**

Storage & Shipping **Store product at  $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.**

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**