

## Anti-AURORA B

Rabbit Polyclonal Antibody

**Catalog # A31-63R**

Lot # B3216-21

### Cited Applications

ELISA, IF, WB

*Ideal working dilutions for each application should be empirically determined by the investigator.*

### Specificity

Recognizes the human Aurora Kinase B protein

### Cross Reactivity

Human

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

The antibody was produced against synthesized peptide corresponding to amino acids 277-238 of Human Aurora Kinase B protein.

### Formulation

0.02 M Potassium Phosphate, 0.15 M Sodium Chloride, pH 7.2 + 0.01% (w/v) Sodium Azide

### Stability

1yr at -20°C from date of shipment

### 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 translocation 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:

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

## Anti-AURORA B

Rabbit Polyclonal Antibody

Catalog #	A31-63R
Lot #	B3216-21
Purification	Immunoaffinity chromatography
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

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.**