

## Aurkb Protein

(Catalog No. E20002)

### Description

AURKB Human is a single, non-glycosylated, polypeptide chain containing 364 amino acids (1-344.. AURKB is fused to 20 a.a. His-Tag at N-terminus and purified by proprietary chromatographic techniques. Aurora Kinase B (AURKB) belongs to a family of mitotic serine/threonine kinases. AURKB connects with chromosomes for the period of prophase prior to relocating to the spindle at anaphase. AURKB localizes to microtubules near kinetochores, specifically to the specialized microtubules called K-fibers. AURKB controls chromosome segregation through the control of microtubule-kinetochore attachment and cytokinesis. AURKB is required for kinetochore localization of BUB1 and SGOL1. AURKB expression during the G2/M phase transition is firmly coordinated with histone H3 phosphorylation, while overexpression is seen in many kinds of cancers. AURKB phosphorylates 'Ser-10' and 'Ser-28' of histone H3 during mitosis. AURKB is a component of the CPC (chromosomal passenger complex), which is a complex that acts as a key regulator of mitosis. High level expression of AURKB is seen in the thymus, which is also expressed in the spleen, lung, testis, colon, placenta and fetal liver. AURKB is expressed during S and G2/M phase and expression is up-regulated in cancer cells during M phase.

### Molecular Weight (kDa)

41.4 kDa

### Formulation

The AURKB solution containing 20mM Tris-HCl buffer (pH8.0), 0.5mM DTT, 20% glycerol, 0.1mM EDTA, 0.1mM EGTA, 0.1M NaCl and 0.1mM PMSF

### Source

Expressed in E.coli expression system

### Size

10 µg

### Storage Temp & Conditions

Store below -18°C. Please prevent freeze thaw cycles. Best use within three months from the date of receipt of this protein.

### Purification Notes

Purity: Greater than 90.0% as determined by SDS-PAGE.

### Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot (Recombinant protein); Antibody Production; Protein Array.

### Ordering Information:

Products	Size	Cat. No.
Aurkb Protein	10 µg	E20002 -1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*