

Aurka Protein

(Catalog No. E20001)

Description

AURKA Human is a single, non-glycosylated, polypeptide chain containing 423 amino acids (1-403). AURKA is fused to 20 a.a. His-Tag at N-terminus and purified by proprietary chromatographic techniques. AURKA (Aurora Kinase A) belongs to the mitotic serine/threonine kinases family. AURKA is a cell cycle-regulated kinase which may be involved in microtubule formation and/or stabilization at the spindle pole during chromosome segregation. AURKA is found at the centrosome in interphase cells and at the spindle poles in mitosis. Since the AURKA expression is cell-cycle regulated, it is low in G1/S, it accumulates during G2/M, and it decreases rapidly after. AURKA is involved in important processes during mitosis and meiosis whose proper function is essential for healthy cell proliferation. In addition, AURKA plays an essential role in tumourigenesis and is overexpressed in various types of cancers. AURKA is strongly expressed in the testis, colon, ovarian, prostate, neuroblastoma, breast and cervical cancer cell lines and weakly in skeletal muscle, thymus and spleen. AURKA interacts with its substrates BORA and ARHGEF2, as well as with TACC1 and CPEB1. Defects in the AURKA gene cause numerical centrosome aberrations including aneuploidy. AURKA overexpression has been linked to chromosomal instability in colorectal cancer. AURKA expression may have a prognostic significance in ovarian carcinoma.

Molecular Weight (kDa)

47.9kDa

Formulation

The AURKA solution containing 20mM Tris-HCl buffer (pH 8.0), 0.5mM DTT, 100mM NaCl, 0.1mM EDTA, 0.1mM EGTA, 0.1mM PMSF and 20% glycerol.

Source

Expressed in E.coli expression system

Size

10 µg

Storage Temp & Conditions

Store below -18°C. Please prevent freeze thaw cycles

Purification Notes

Purity: Greater than 95.0% as determined by SDS-PAGE

Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot (Recombinant protein); Antibody Production; Protein Array; Preparation Method: in vitro wheat germ expression system ;Best use within three months from the date of receipt of this protein

Ordering Information:

Products	Size	Cat. No.
Aurka Protein	10 µg	E20001-1

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