

AURORA C, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog # A33-14G

Lot # X575-1

Product Description

Recombinant full-length human AURORA C was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_003160](#). It can be activated by INCENP *in vitro*.

Gene Aliases

AURKC, AIE2, AIK3, AurC, STK13, aurora-C

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°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.

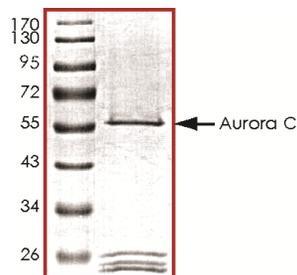
Scientific Background

AURORA C is a member of mitotic serine/threonine kinases that regulate centrosome maturation, chromosome segregation, and cytokinesis. AURORA C has been shown to contribute to oncogenic transformation and is frequently overexpressed and amplified in many human tumors (1). AURORA C overexpression has been correlated with centrosome amplification, which can be a driving cause of genomic instability in tumor cells. AURORA C is an important regulator of both genomic integrity and cell cycle progression in cancer cells and is an attractive target for anticancer drug development (2).

References

1. Steven, L. et al: Targeting Aurora-2 Kinase in Cancer. *Molecular Cancer Therapeutics*. 2003;2: 589-595.
2. Shingo, T. et al: The centrosomal protein Lats2 is a phosphorylation target of Aurora-A kinase. *Genes to Cells*. 2004; 9: 383-397.

Purity



The purity of Aurora C was determined to be **>70%** by densitometry. Approx. MW **61kDa**.

AURORA C, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	A33-14G
Specific Lot Number	X575-1
Purity	>70%
Concentration	0.1 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°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
www.signalchem.com

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