

Anti-BUB1

Rabbit Polyclonal Antibody

Catalog # B11-63R

Lot # O2121-6

Cited Applications

WB, ELISA, ICC, IF

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

Specificity

Recognizes the BUB1 protein

Cross Reactivity

Human and Mouse

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

BUB1 antibody was raised against a 19 amino acid synthetic peptide from near the amino terminus of human BUB1

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

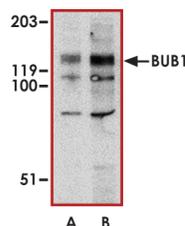
Scientific Background

BUB1 is a kinase which involved in spindle checkpoint function that in part by phosphorylating a member of the mitotic checkpoint complex and activating the spindle checkpoint. BUB1 is highly expressed in thymus, ovary, and testis (1). BUB1 maintains the steady-state level and centromeric localization of SGO1, and they concluded that BUB1 protects centromeric cohesion in mitosis through SGO1 (2). BUB1-mediated phosphorylation of CDC20 is required for proper checkpoint signaling and that inhibition of APC by BUB1 may partly account for the sensitivity of the spindle checkpoint. BUB1 kinase creates a mark for shugoshin localization and the correct partitioning of chromosomes.

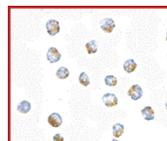
References

1. Leland, S.et.al: Heterozygosity for a Bub1 mutation causes female-specific germ cell aneuploidy in mice. Proc. Nat. Acad. Sci. 106: 12776-12781, 2009.
2. Tang, Z.et.al: Human Bub1 protects centromeric sister-chromatid cohesion through Shugoshin during mitosis. Proc. Nat. Acad. Sci. 101: 18012-18017, 2004.

Sample Data



Western blot analysis of BUB1 in A-20 lysate with BUB1 antibody at (A) 2 and (B) 4 ug/ml.



Immunocytochemistry of BUB1 in A20 cells with BUB1 antibody at 10 ug/ml.

Anti-BUB1

Rabbit Polyclonal Antibody

Catalog Number

B11-63R

Specific Lot Number

O2121-6

Purification

Affinity 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
www.signalchem.com

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