

Anti-BUB1B

Mouse Monoclonal Antibody

Catalog # B12-60M

Lot # B2316-24

Cited Applications

ELISA, IF, IP, WB

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

Specificity

Recognizes the human BUBR1 Kinase protein

Cross Reactivity

Human

Host/Isotype/Clone#

Mouse, IgG1

Immunogen

The antibody was produced against synthesized peptide corresponding to amino acid residues 1-350 of human BUBR1 Kinase 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

BUB1B or budding uninhibited by benzimidazoles 1 homolog beta is involved in spindle checkpoint function and has been localized to the kinetochore. BUB1B plays a role in the inhibition of the anaphase-promoting complex/cyclosome (APC/C), delaying the onset of anaphase and ensuring proper chromosome segregation (1). BUB1B links the regulation of chromosome-spindle attachment to mitotic checkpoint signaling (2). BUB1B expression during the spindle checkpoint is dependent on residual BUB1B expression and impaired spindle checkpoint function has been found in many forms of cancer. BUB1B is essential for early embryonic development and normal hematopoiesis.

References:

1. Rio Frio. Et.al: Homozygous BUB1B mutation and susceptibility to gastrointestinal neoplasia. *New Eng. J. Med.* 363: 2628-2637, 2010.
2. Lampson, M. A., Kapoor. et.al: The human mitotic checkpoint protein BubR1 regulates chromosome-spindle attachments. *Nature Cell Biol.* 7: 93-98, 2005.

Anti-BUB1B

Mouse Monoclonal Antibody

Catalog #	B12-60M
Lot #	B2316-24
Purification	Protein A 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.