

Bub3 (CT) antibody (pAb)

Rabbit Anti-Human/Mouse Bub3 (CT)

Instruction Manual

Catalog Number	PK-AB718-4227
Synonyms	Bub3 Antibody: Budding uninhibited by benzimidazoles 3, Bub3L
Description	The mitotic checkpoint protein Bub3 is involved with the essential spindle checkpoint pathway which operates during early embryogenesis. Bub3 is important during G2 and early mitosis stages, permitting entry into mitosis depending upon the assembly state of microtubules, thus preventing premature sister chromatid separation, mis-segregation and aneuploidy. Bub3 contains four WD repeat domains and is required for the kinetochore localization of Bub1, a related kinase that is necessary for spindle assembly checkpoint function. Bub1 is able to autophosphorylate and can catalyze the phosphorylation of Bub3. Both Bub1 and Bub3 are mutually dependent for function. Altered Bub expression levels may significantly impair mitotic checkpoint function and is associated with tumor cell proliferation.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Bub3 antibody was raised in rabbits against a 16 amino acid peptide from near the carboxy terminus of human bub3.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, ICC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Bub3 antibody can be used for detection of bub3 by Western blot at 0.5 - 1 µg/mL. Antibody can also be used for immunocytochemistry starting at 10 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Kalitsis P, Earle E, Fowler KJ, et al. Bub3 gene disruption in mice reveals essential mitotic spindle checkpoint function during early embryogenesis. <i>Genes Dev.</i> 2000; 14:2277-82. Taylor SS, Ha E and McKeon F. The human homologue of Bub3 is required for kinetochore localization of Bub1 and a Mad3/Bub1-related protein kinase. <i>J. Cell Biol.</i> 1998; 142:1-11. Warren CD, Brady DM, Johnston RC, et al. Distinct chromosome segregation roles for spindle checkpoint proteins. <i>Mol. Biol. Cell</i> 2002; 13: 3029-41. Roberts BT, Farr KA and Hoyt MA. The <i>Saccharomyces cerevisiae</i> checkpoint gene BUB1 encodes a novel protein kinase. <i>Mol. Cell Biol.</i> 1994; 14:8282-91.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4227P; Bub3 Peptide Cat.No. PK-AB718-1205; Jurkat Cell Lysate

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.