

Tubulin antibody (pAb)

Rabbit Anti-Human/Mouse/Rat/Dog Tubulin

Instruction Manual

Catalog Number	PK-AB577-3708
Description	Microtubules are one type of cytoskeleton. Globular tubulin subunits are the building blocks of a microtubule and the alpha and beta-tubulin heterodimer forms the tubulin subunits, which is present in all eukaryotic cells. Microtubules mediate many cell movements.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide corresponding to the C-terminus of human beta-Tubulin.
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human, mouse, rat, dog
Specificity	The antibody recognizes 55 kDa beta-tubulin in samples of human, mouse, and rat origins. Reactivity to other species has not been determined.
Formulation	100 µg (0.5 mg/ml) affinity purified rabbit polyclonal antibody in phosphate-buffered saline (PBS) containing 30% glycerol, 0.5% BSA, and 0.01% thimerosal.
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Storage & Stability	Store at -20°C. For long-term storage, aliquot and refreeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	Western blotting (0.5-4 µg/ml). The optimal concentrations should be determined individually. Antibody might be suitable for other applications not tested so far. HeLa cell or Jurkat cell lysate can be used as a positive control.

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