

HEC-1 antibody (pAb)

Rabbit Anti-Human/Mouse HEC-1

Instruction Manual

Catalog Number	PK-AB577-3763
Description	HEC, HEC1, HEC-1, Highly Expressed in Cancer 1, KNTC2, Kinetochore Associated 2
Quantity	100 µl
Source / Host	Rabbit
Immunogen	Synthetic peptide of human HEC1
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human, mouse
Specificity	See Applications or Species Reactivity.
Formulation	100 µl affinity purified rabbit polyclonal antibody in phosphate-buffered saline (PBS) containing 30% glycerol, 1% BSA and 0.02% 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 (1:500 – 1:2500) and immunoprecipitation. HeLa and NIH3T3 cell lysates can be used as a positive control. The optimal concentrations should be determined individually. Antibody might be suitable for other applications not tested so far. The antibody recognizes the HEC1 of human and mouse origins. Reactivity to other species has not been tested.

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