

Instruction Manual

Catalog Number	PK-AB577-3707
Description	Histone acetyltransferases (HATs) have been implicated in a number of cellular functions including gene regulation, DNA synthesis, and repair. Histone acetyltransferases and deacetylases are respectively, the enzymes devoted to the addition and removal of acetyl groups from lysine residues on the Histone N-terminal tails. The enzymes exert fundamental roles in developmental processes and their deregulation has been linked to the progression of diverse human disorders, including cancer.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide containing amino acids near the C-terminus of the human HAT-3 (ID BV-85).
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	The antibody recognizes 220 kDa human HAT-3. A 50 kDa unknown band can also be detected. Reactivity to other species has not been determined.
Formulation	100 µg (0.5 mg/ml) Protein A 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 freeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	The antibody can be used for Western blotting (2-6 µg/ml). However, the optimal concentrations should be determined individually. Jurkat cell lysate (Cat.# PK-AB718-1205) can be used as a positive control. Blocking peptide is available separately (Cat.# PK-AB577-3707P).

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