

Instruction Manual

Catalog Number	PK-AB577-3689
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 surrounding amino acid 445 of human HAT-1 (ID BV-249).
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	The antibody recognizes 50 kDa HAT-1 from samples of human, mouse and rat origins.
Formulation	100 µg (0.2 mg/ml) immunoaffinity 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 (1-2 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 (Cat.# PK-AB577-3689P) is also available.

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