

Instruction Manual

Catalog Number	PK-AB577-3605
Description	Human HDAC-5 is composed of 1122 amino acid residues. The deacetylase domain of HDAC-5 is located at the C-terminal half of the molecule. The N-terminal non-deacetylase domain does not show any significant homology with any published sequence. Both domains are required for HDAC-5-mediated repression of gene transcription. HDAC-5 interacts with a growing number of transcriptional factors including MEF2A as well as other HDAC proteins. The interacting complexes bind to specific regions of chromatin and regulate gene transcription in these regions.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide surrounding amino acid 575 of human HDAC-5 (Internal ID BV-78).
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	The antibody detects ~124 kDa histone deacetylase 5. It does not cross-react with other HDAC proteins including HDAC-1, 2, 3, 4, 6, 7, 8, 9,10, and 11.
Formulation	100 µg (0.5 mg/ml) affinity purified rabbit anti-HDAC-5 polyclonal antibody in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 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 (0.5-4 µg/ml). Based on researchers feedback, the antibody also works well in Immunoprecipitation (20 µg/ml), and Immunohistochemistry (20 µg/ml). However, the optimal conditions should be determined individually. Blocking peptide (Cat.# PK-AB577-3605P) is available separately.

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