

## Instruction Manual

Catalog Number	PK-AB577-3603
Description	HDAC-3 is a member of the class I mammalian histone deacetylase family, which plays an important role in modulating the eukaryotic chromatin structure. Human HDAC-3 is composed of 428 amino acid residues. HDAC-3 interacts with a growing number of transcriptional factors including SMRT and N-CoR. 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 corresponding to the C-terminus of human HDAC-3 (ID BV-224).
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat, Chicken
Specificity	The antibody detects the 49 kDa HDAC-3. It does not cross-react with other HDAC proteins including HDAC-1, 2, 4, 5, 6, 7, and 8.
Formulation	100 µg (0.2 mg/ml) peptide affinity purified rabbit anti-HDAC-3 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, it can also be used in Immunohistochemistry (20-30 µg/ml). However, the optimal conditions should be determined individually. Jurkat cell lysate (Cat.# PK-AB718-1205) can be used as a positive control. Blocking peptide (Cat.# PK-AB577-3603P) is also available.

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