

Instruction Manual

Catalog Number	PK-AB577-3601
Description	HDAC-1 is an active component of transcriptional corepressor complexes which can be recruited to specific promoter regions via transcriptional factors. HDAC-1 catalyzes removal of acetyl-groups from acetyl-lysines of histones and promotes compaction of chromatin in these regions, which leads to the inhibition of gene transcription.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide corresponding to the C-terminus of human HDAC1 (ID BV-220).
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	The antibody detects the 62 kDa histone deacetylase 1. It does not cross-react with other HDAC proteins including HDAC-2, 3, 4, 5, 6, 7, and 8.
Formulation	100 µg (0.5 mg/ml) peptide affinity purified rabbit anti-HDAC-1 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 researcher's feedback, it also works well in Immunoprecipitation (10-20 µg/ml), and Immunohistochemistry (10-20 µg/ml). However, the optimal conditions should be determined individually. Blocking peptide is available separately (Cat.# PK-AB577-3601P).

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