

Instruction Manual

Catalog Number	PK-AB577-3610
Description	HDAC family are divided into two classes, I and II. Class I of the HDAC family comprises four members, HDAC-1, 2, 3, and 8. Class II of the HDAC family comprises HDAC-4, 5, 6, and 7, the molecular weights of which are all about two-fold larger than those of the class I members. Human HDAC-1, 2 and 3 were expressed in various tissues, but the others (HDAC-4, 5, 6, and 7) showed tissue-specific expression patterns. These results suggest that each member of the HDAC family exhibits a different, individual substrate specificity and function in vivo.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide near the N-terminus of human HDAC-10 (ID BV-73).
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	The antibody recognizes 74 kDa HDAC-10 of human, mouse, and rat origins. A 55 kDa cleavage fragment can also be detected in mouse and rat tissue lysates.
Formulation	100 µg (0.2 mg/ml) 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 in Western blotting (0.5-4 µg/ml), However, the optimal concentrations should be determined individually. Blocking peptide (Cat.# PK-AB577-3610P) is available separately.

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