

HDAC-2 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat/Chicken/Hamster/Dog Histone Deacetylase 2

Instruction Manual

Catalog Number	PK-AB577-3602
Description	HDAC-2 is a class I mammalian histone deacetylase containing 488 amino acid residues with an apparent size of ~55 kDa by denaturing SDS-PAGE. HDAC-2 has been shown to interact directly with transcription repressor complexes and the nuclear receptor corepressor N-CoR.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide surrounding amino acid 479 of human HDAC2 (ID BV-222).
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat, Chicken, Hamster, Dog
Specificity	The antibody detects 55 kDa Histone Deacetylase 2 in samples from human, mouse, rat, chicken, hamster, and canine origins.
Formulation	100 µg (0.2 mg/ml) affinity purified rabbit anti-HDAC-2 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 Immunocytochemistry (10-20 µg/ml). However, the optimal conditions should be determined individually. Blocking peptide is available separately (Cat.# PK-AB577-3602P).

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