

HDAC-8 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Histone Deacetylase 8

Instruction Manual

Catalog Number	PK-AB577-3608
Description	Mammalian HDAC-8, isolated from human kidney, is a member of the histone deacetylase family. HDAC-8 shares homology with other HDACs in the family, but has different tissue distribution. HDAC-8 is localized to the nucleus and plays a role in the tissue development as well as cancer etiology.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide surrounding amino acid 10 of mouse HDAC-8 (ID BV-70).
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	The antibody recognizes 78 kDa and 42 kDa HDAC-8 of human, mouse, and rat origins.
Formulation	100 µg (0.2 mg/ml) protein A 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 has been used in Western blotting (0.5-4 µg/ml). Based on researchers feedback, it also works well in immunoprecipitation (10-20 µg/ml), and immunohistochemistry (10-20 µg/ml). However, the optimal concentrations should be determined individually. Blocking peptide (Cat.# PK-AB577-3608P) is also available separately.

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