

HDAC-7 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Histone Deacetylase 7

Instruction Manual

Catalog Number	PK-AB577-3607
Synonyms	HDAC7A, histone deacetylase 7A
Description	Human HDAC-7 is composed of 912 amino acid residues. Although HDAC-7 is localized mostly to the cell nucleus, it is also found in the cytoplasm, suggesting nucleo-cytoplasmic shuttling. The histone deacetylase activity of HDAC-7 maps to a carboxy-terminal domain and is dependent on interaction with class I HDACs in the nucleus. HDAC-7 catalyzes removal of acetyl-groups from acetyl-lysines of histones and promotes compaction of chromatin in these regions, leading to the inhibition of gene transcription.
Quantity	100 µg
Source / Host	Rabbit
Subclass	Rabbit IgG
Immunogen	Synthetic peptide surrounding amino acid 921 of murine HDAC7 (ID BV-79).
Specific Species React.	Human, Mouse, Rat
Cross-Reactivity	The antibody detects ~105 kDa HDAC-7.
Formulation	100 µg (0.5 mg/ml) affinity purified rabbit anti-HDAC-7 polyclonal antibody in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 5 mM EDTA, 0.03% Proclin.
Stabilizer	thimerosal
Concentration	0.5 mg/ml
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Applications	The antibody can be used in Western blotting (0.5-4 µg/ml). Based on researchers feedback, it can also be used in immunoprecipitation (20-40 µg/ml) and immunohistochemistry (20-40 µg/ml). However, the optimal conditions should be determined individually. Mouse small intestine tissue lysate (Cat.# PK-AB718-1408) can be use as a positive control. Blocking peptide (Cat.# PK-AB577-3607P) is also available.
Storage	Store at -20°C. For long-term storage, aliquot and freeze at -70°C. Avoid repeated freeze/thaw cycles.
Shipping Conditions	2-8°C

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