

# HDAC-4 antibody (pAb)

## Rabbit Anti-Human/Mouse/Rat Histone Deacetylase 4

### Instruction Manual

Catalog Number	PK-AB577-3604
Description	HDAC-4 is a member of the class II mammalian histone deacetylases, which consists of 1084 amino acid residues. Its C-terminal sequence is highly similar to the deacetylase domain of yeast HDA-1. HDAC-4, unlike other deacetylases, shuttles between the nucleus and cytoplasm in a process involving active nuclear export. Association of HDAC-4 with 14-3-3 results in sequestration of HDAC-4 protein in the cytoplasm. In the nucleus, HDAC-4 associates with the myocyte enhancer factor MEF2A. Binding of HDAC-4 to MEF2A results in the repression of MEF2A transcriptional activation. HDAC-4 has also been shown to interact with other deacetylases such as HDAC-3 as well as the corepressors NcoR and SMART.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	NA
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	The antibody detects ~140 kDa HDAC-4. It does not cross-react with other HDAC proteins including HDAC-1, 2, 3, 5, 6, 7, and 8.
Formulation	100 µg (0.5 mg/ml) affinity purified rabbit anti-HDAC-4 polyclonal antibody in phosphate buffered saline (PBS), pH 7.2, containing 50% glycerol, 1% BSA, 0.02% sodium azide.
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 (1 µg/ml). However, the optimal conditions should be determined individually. Blocking peptide (Cat.# PK-AB577-3604P) is available separately.

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