

# HDAC-11 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Histone Deacetylase 11

## Instruction Manual

<b>Catalog Number</b>	PK-AB577-3611
<b>Synonyms</b>	HD11, Histone deacetylase 11
<b>Description</b>	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.
<b>Quantity</b>	100 µg
<b>Source / Host</b>	Rabbit
<b>Subclass</b>	Rabbit IgG
<b>Immunogen</b>	Synthetic peptide surrounding amino acid 269 of human HDAC-11 (ID BV-75).
<b>Specific Species React.</b>	Human, Mouse, Rat
<b>Cross-Reactivity</b>	The antibody recognizes 39 kDa HDAC-11 of human, mouse, and rat origins. APPLICATION: The antibody can be used in Western blotting (0.5-4 µg/ml), However, the optimal concentrations should be determined individually.
<b>Formulation</b>	0.5 mg/ml in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 5 mM EDTA and 0.035 Proclin.
<b>Reconstitution</b>	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
<b>Applications</b>	The antibody can be used in Western blotting (0.5-4 µg/ml), However, the optimal concentrations should be determined individually. Blocking peptide is available separately (Cat.# PK-AB577-3611P).
<b>Storage</b>	Store at -20°C. For long-term storage, aliquot and freeze at -70°C. Avoid repeated freeze/thaw cycles.
<b>Shipping Conditions</b>	2-8°C

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