

## HDAC6 Polyclonal Antibody

(Catalog # A-4006)

### Background

Histone deacetylases (HDACs) play a critical role in transcriptional repression of gene expression in eukaryotic cells through catalyzing the hydrolytic removal of acetyl groups from histone lysine residues. Three distinct families of HDACs have been described, comprising a group of at least 20 proteins in humans. HDAC6 is a class II histone deacetylase.

### Description

HDAC6 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

### Specificity

Human, Mouse, Rat

### Isotype

IgG

### Uniprot ID

Q9UBN7

### Purification

Affinity Purified

### Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-200 of human HDAC6 (NP\_006035.2).

### Storage

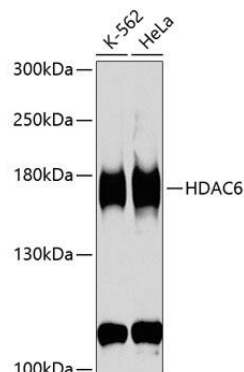
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

### Alternative Names

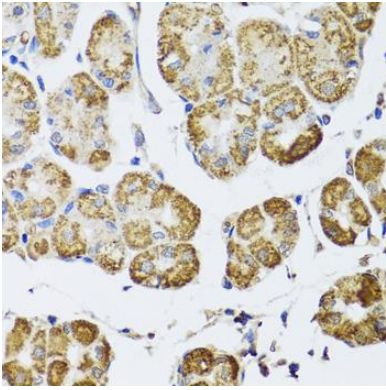
FLJ16239 antibody, HD 6 antibody, Histone deacetylase 6 (HD6) antibody, JM 21 antibody, KIAA0901 antibody, OTTHUMP00000032398 antibody, OTTHUMP00000197663 antibody

### Application

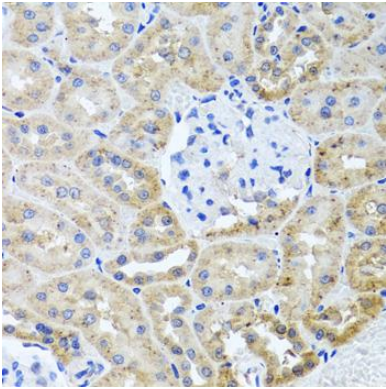
WB, IHC; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:200



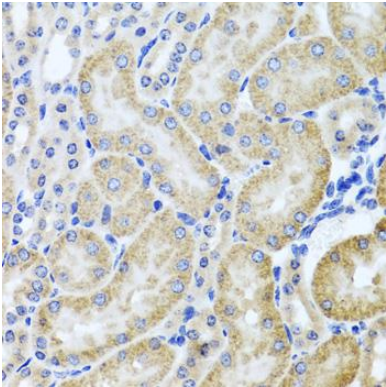
Western blot analysis of extracts of various cell lines, using HDAC6 antibody at 1:1000 dilution.  
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.  
Lysates/proteins: 25ug per lane.  
Blocking buffer: 3% nonfat dry milk in TBST.



Immunohistochemistry of paraffin-embedded human stomach using HDAC6 antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded rat kidney using HDAC6 antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded mouse kidney using HDAC6 antibody at dilution of 1:100 (40x lens).