

---

## HDAC4 Polyclonal Antibody

(Catalog # A70941)

---

### Background

Histones play a critical role in transcriptional regulation, cell cycle progression, and developmental events. Histone acetylation/deacetylation alters chromosome structure and affects transcription factor access to DNA. The protein encoded by this gene belongs to class II of the histone deacetylase/acuc/apha family. It possesses histone deacetylase activity and represses transcription when tethered to a promoter. This protein does not bind DNA directly, but through transcription factors MEF2C and MEF2D. It seems to interact in a multiprotein complex with RbAp48 and HDAC3.

### Description

HDAC4 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Human, Mouse, Rat

### Isotype

IgG

### Uniprot ID

P56524

### Purification

Affinity Purification

### Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 451-650 of human HDAC4 (NP\_006028.2).

### Storage

Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

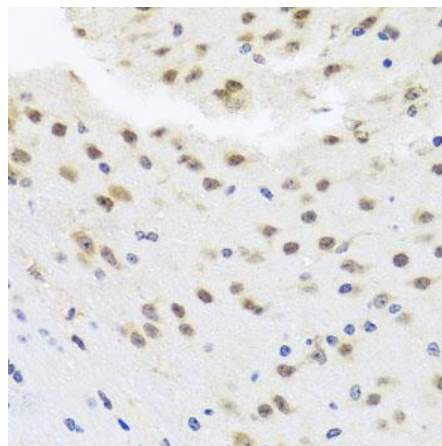
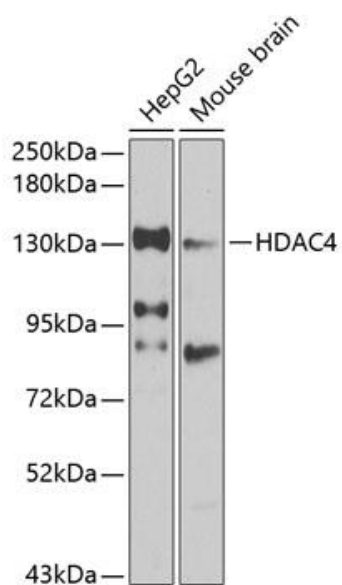
### Alternative Names

HDAC4; AHO3; BDMR; HA6116; HD4; HDAC-4; HDAC-A; HDACA; histone deacetylase 4

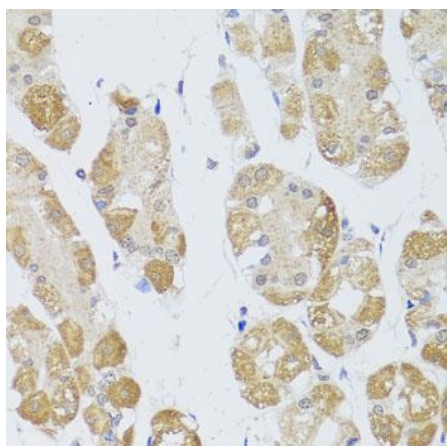
### Application

WB, IHC; Recommended dilution: WB 1:500 - 1:1000, IHC 1:50 - 1:100

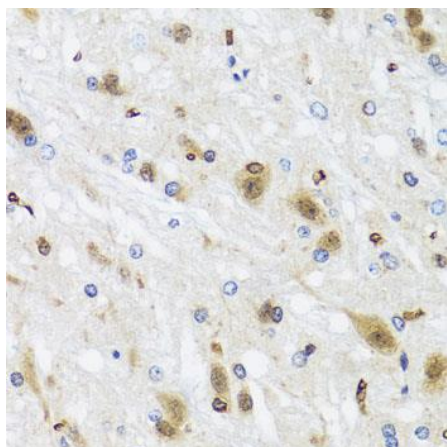
---



Immunohistochemistry of paraffin-embedded mouse brain using HDAC4 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded human stomach using HDAC4 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded rat brain using HDAC4 Polyclonal Antibody at dilution of 1:100 (40x lens).