

## HDAC10 Polyclonal Antibody

(Catalog # A-4010)

### 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. HDAC10 is a novel class II histone deacetylase.

### Description

HDAC10 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

0.03% Proclin 300. 50% Glycerol, 0.01M PBS, PH 7.4.

### Specificity

Human, Mouse

### Isotype

IgG

### Uniprot ID

Q969S8

### Purification

Protein G purified

### Immunogen

Recombinant Human Histone deacetylase 10 protein (207-412AA)

### Storage

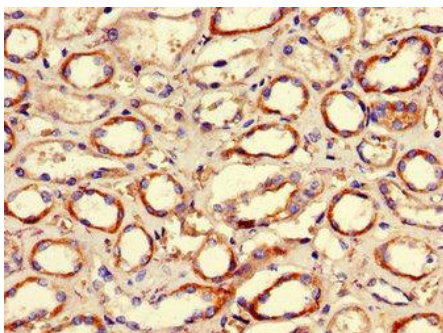
Shipped at 4°C (Blue Ice). Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

### Alternative Names

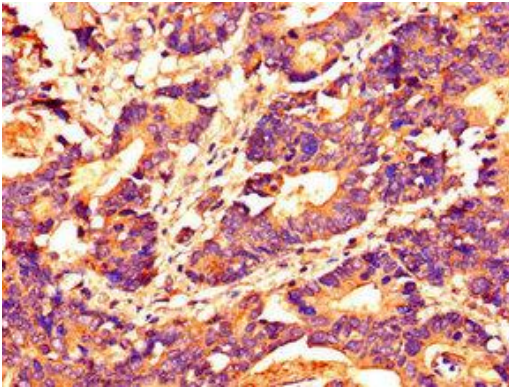
DKFZP761B039 antibody, HD 10 antibody, Histone deacetylase 10 antibody, MGC149722 antibody, OTTHUMP00000028555 antibody

### Application

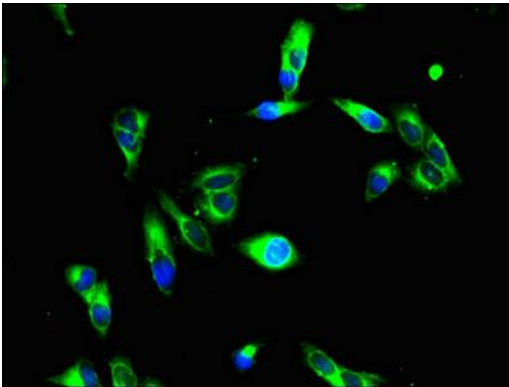
ELISA, WB, IHC, IF; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200, IF:1:50-1:200



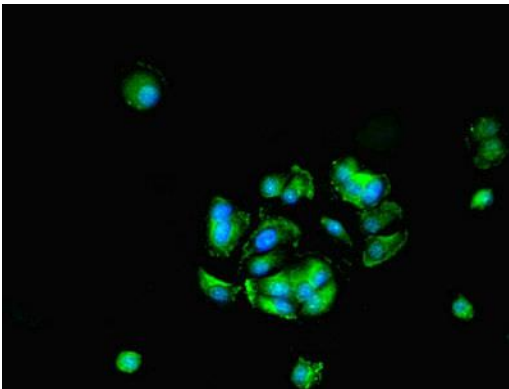
Immunohistochemistry of paraffin-embedded human kidney tissue at dilution of 1:100



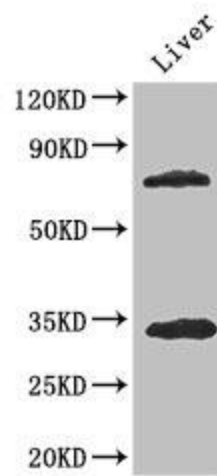
Immunohistochemistry of paraffin-embedded human colon cancer at dilution of 1:100



Immunofluorescent analysis of HeLa cells at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunofluorescent analysis of MCF-7 cells at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Western Blot  
Positive WB detected in: Mouse liver tissue  
All lanes: HDAC10 Polyclonal antibody at 3.4µg/ml  
Secondary  
Goat polyclonal to rabbit IgG at 1/50000 dilution  
Predicted band size: 72, 70, 71, 44 kDa  
Observed band size: 72, 34 kDa