

HDAC9 Recombinant Monoclonal Antibody [1F2]

(Catalog # A73829)

Background

Responsible for the deacetylation of lysine residues on the N-terminal part of the core histones (H2A, H2B, H3 and H4). Histone deacetylation gives a tag for epigenetic repression and plays an important role in transcriptional regulation, cell cycle progression and developmental events. Represses MEF2-dependent transcription.

Description

HDAC9 Recombinant Monoclonal Antibody [1F2]. Unconjugated. Raised in: HEK293F Cell.

Formulation

Buffer: Rabbit IgG in phosphate buffered saline, pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

Specificity

Human

Isotype

Rabbit IgG

Uniprot ID

Q9UKV0

Purification

Affinity Chromatography

Immunogen

A synthesized peptide derived from human HDAC9

Storage

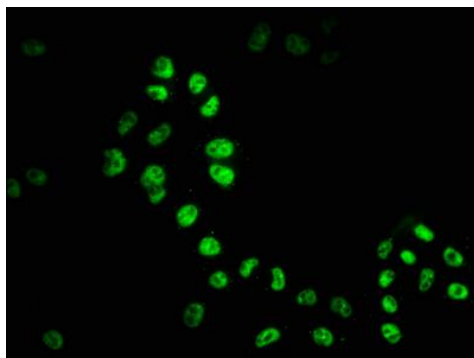
Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

Alternative Names

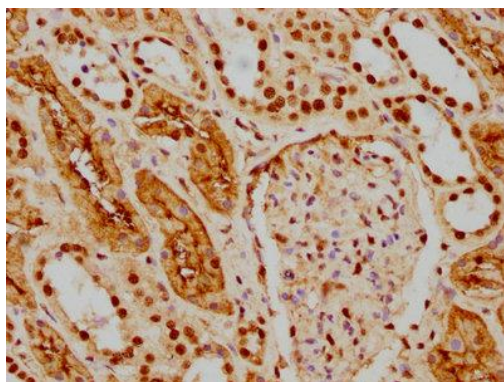
Histone deacetylase 9, HD9, Histone deacetylase 7B, HD7, HD7b, Histone deacetylase-related protein, MEF2-interacting transcription repressor MITR, HDAC9, HDAC7, HDAC7B, HDRP, KIAA0744, MITR

Application

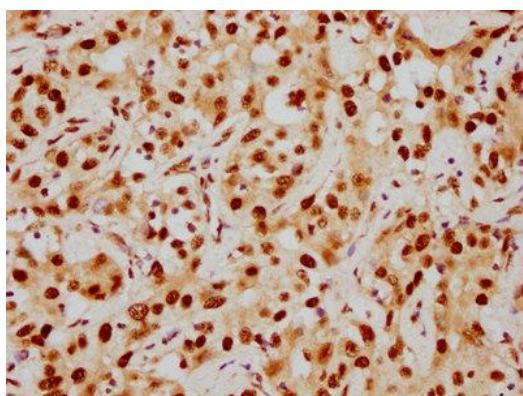
ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:50-1:200, IF:1:20-1:200



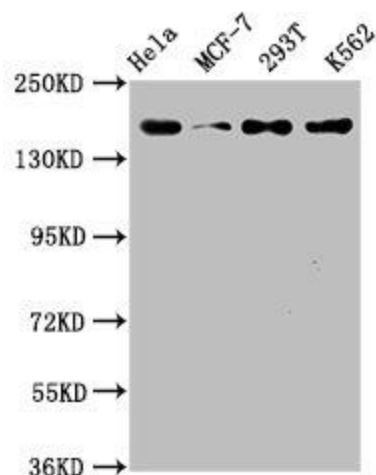
Immunofluorescence staining of HeLa cells with HDAC9 rMAb at 1:51, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L).



IHC image of HDAC9 rMAb diluted at 1:154 and staining in paraffin-embedded human kidney tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



IHC image of HDAC9 rMAb diluted at 1:154 and staining in paraffin-embedded human liver cancer performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



Western Blot

Positive WB detected in: HeLa whole cell lysate, MCF-7 whole cell lysate, 293T whole cell lysate, K562 whole cell lysate

All lanes: HDAC9 antibody at 1.54ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 112, 102, 66, 98, 118, 113, 61, 63, 58 KDa. Observed band size: 160 KDa