
HDAC3 Polyclonal Antibody

(Catalog # A68305)

Background

Responsible for the deacetylation of lysine residues on the N-terminal part of the core histones (H2A, H2B, H3 and H4), and some other non-histone substrates. Histone deacetylation gives a tag for epigenetic repression and plays an important role in transcriptional regulation, cell cycle progression and developmental events. Histone deacetylases act via the formation of large multiprotein complexes. Participates in the BCL6 transcriptional repressor activity by deacetylating the H3 'Lys-27' (H3K27) on enhancer elements, antagonizing EP300 acetyltransferase activity and repressing proximal gene expression. Probably participates in the regulation of transcription through its binding to the zinc-finger transcription factor YY1; increases YY1 repression activity. Required to repress transcription of the POU1F1 transcription factor. Acts as a molecular chaperone for shuttling phosphorylated NR2C1 to PML bodies for sumoylation.

Description

HDAC3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

O15379

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Histone deacetylase 3 protein (1-428AA)

Storage

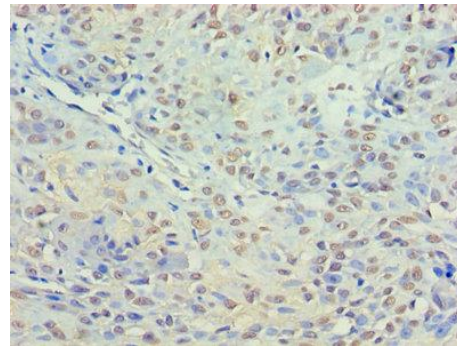
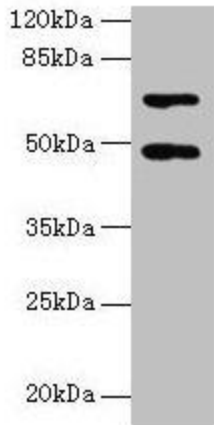
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

HD3, RPD3-2, SMAP45, HDAC3

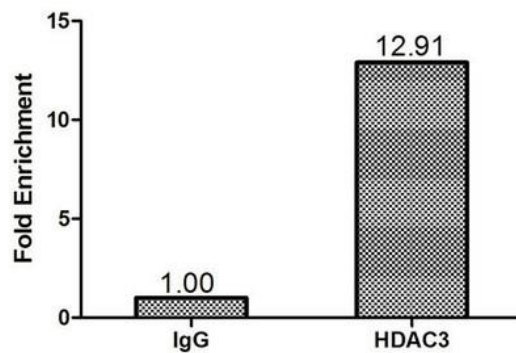
Application

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

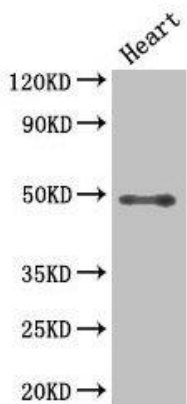


Immunohistochemistry of paraffin-embedded human breast cancer using HDAC3 Antibody at dilution of 1:100

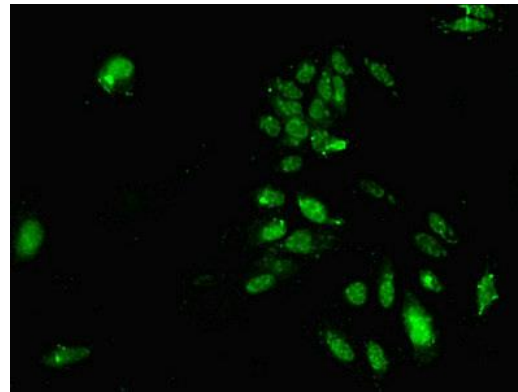
Western blot
 All lanes: HDAC3 antibody at 2ug/ml + HeLa whole cell lysate
 Secondary
 Goat polyclonal to rabbit IgG at 1/10000 dilution
 Predicted band size: 49, 50 kDa
 Observed band size: 49 kDa



Chromatin Immunoprecipitation HeLa (1.2×10^6) were cross-linked with formaldehyde, sonicated, and immunoprecipitated with 4ug anti-HDAC3 or a control normal rabbit IgG. The resulting ChIP DNA was quantified tissue using real-time PCR with primers against the P21 promoter.



Western Blot
 Positive WB detected in: Mouse heart tissue
 All lanes: HDAC3 antibody at 2.5ug/ml
 Secondary
 Goat polyclonal to rabbit IgG at 1/50000 dilution
 Predicted band size: 49, 50 kDa
 Observed band size: 49 kDa



Immunofluorescent analysis of HeLa cells using HDAC3 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)