
HDAC5 Polyclonal Antibody

(Catalog # A-4005)

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. HDAC5 is a class II histone deacetylase.

Description

HDAC5 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse, Rat, Human

Isotype

IgG

Uniprot ID

Q9UQL6

Purification

Affinity Purified

Immunogen

Synthetic Peptide of Human HDAC5a-3018

Storage

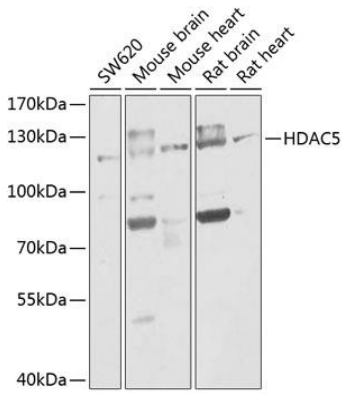
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

Antigen NY CO 9 antibody, FLJ90614 antibody, HD 5 antibody, Histone deacetylase 5 antibody, KIAA0600 antibody

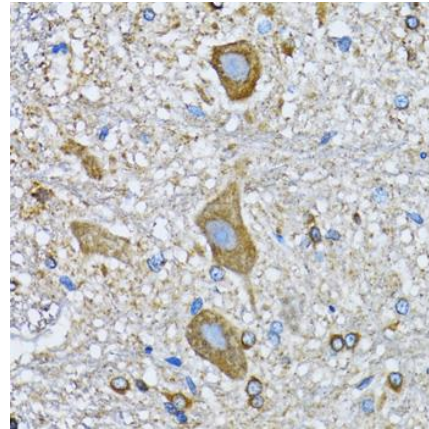
Application

WB, IHC, IF; Recommended dilution: WB 1:500 - 1:2000, IHC-P 1:50 - 1:200, IF/ICC 1:50 - 1:200

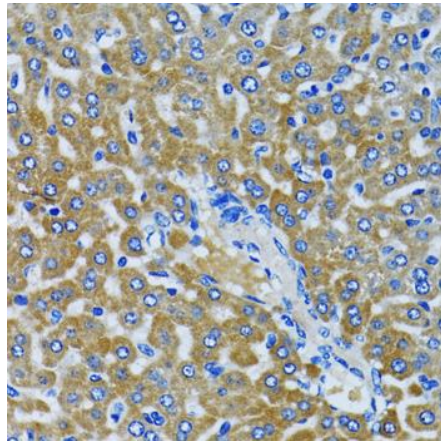
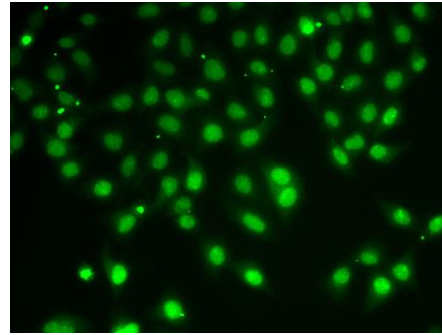


Western blot analysis of extracts of various cell lines, using HDAC5 antibody at 1:1000 dilution.
 Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25ug per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.

Immunohistochemistry of paraffin-embedded human tonsil using HDAC5 antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded mouse spinal cord using HDAC5 antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded rat liver using HDAC5 antibody at dilution of 1:100 (40x lens).

Immunofluorescence analysis of A-549 cells using HDAC5 antibody.

