

HDAC7 Polyclonal Antibody

(Catalog # A-4007)

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. HDAC7 is a class II histone deacetylase.

Description

HDAC7 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

Q8WUI4

Purification

Affinity Purified

Immunogen

Synthetic Peptide of Human HDAC7

Storage

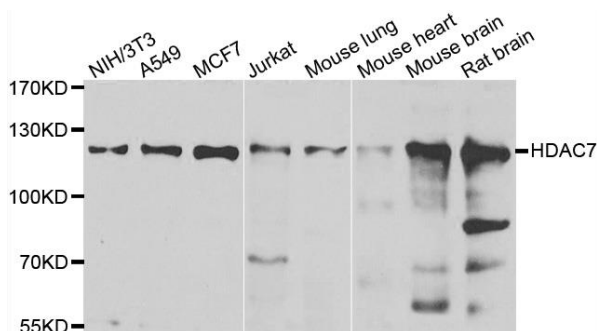
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

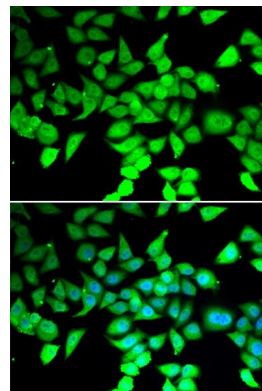
DKFZP586J0917 antibody, FLJ99588 antibody, HD7 antibody, HD7a antibody, HDAC 7A antibody, Histone deacetylase 7 antibody, Histone deacetylase 7A antibody, OTTHUMP00000202813 antibody, OTTHUMP00000202814 antibody

Application

WB, IF



Western blot analysis of extracts of various cell lines, using HDAC7 Polyclonal Antibody.



Immunofluorescence analysis of MCF7 cell using HDAC7 Polyclonal Antibody. Blue: DAPI for nuclear staining.