

HDAC3 Antibody

Cat. No. C15200145

Type: Monoclonal	Specificity: Human
Size: 50 µg/25 µl	Isotype: IgG1
Concentration: 2.0 µg/µl	Host: Mouse
Lot No.: 001	Purity: Protein A purified monoclonal antibody
Storage buffer: PBS containing 0.05% azide and 0.05% ProClin 300.	Storage conditions: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.
Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.	

Description

Alternative names: **HD3, RPD3, RPD3-2, SMAP45**

Monoclonal antibody raised in mouse against human **HDAC3 (Histone deacetylase 3)**, using a KLH-conjugated synthetic peptide containing a sequence from the C-terminal region of the protein.

Applications

Applications	Suggested dilution	References
Immunofluorescence	1:500	Fig 1

Target Description

HDAC3 (UniProt/Swiss-Prot entry O15379) catalyses the deacetylation of lysine residues in the N-terminal part of the core histones (H2A, H2B, H3 and H4). Acetylation and deacetylation of these highly conserved lysine residues is important for the control of gene expression and HDAC activity is associated with gene repression. Histone deacetylation is established by the formation of large multiprotein complexes. HDAC3 may bind to the zinc-finger transcription factor YY1, thereby regulating transcription. It is also able to modulate cell growth and apoptosis through the interaction with p53 and is thought to be essential for the repression of the POU1F1 transcription factor.

Validation Data

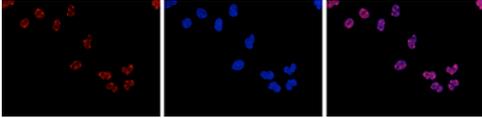


Figure 1. Immunofluorescence using the Diagenode monoclonal antibody directed against HDAC3

HeLa cells were stained with the Diagenode antibody against HDAC3 (Cat. No. C15200145) and with DAPI. Cells were fixed with 4% formaldehyde for 10' and blocked with PBS/TX-100 containing 5% normal goat serum and 1% BSA. The cells were immunofluorescently labelled with the HDAC3 antibody (left) diluted 1:500 in blocking solution followed by an anti-mouse antibody conjugated to Alexa594. The middle panel shows staining of the nuclei with DAPI. A merge of the two stainings is shown on the right.