

HDAC5 polyclonal antibody - Classic

Cat. No. C15410275

Type: Polyclonal

Source: Rabbit

Lot #: 001

Size: 100 µg

Concentration: 1.12 µg/µl

Specificity: Human

Purity: Affinity purified

Storage: 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.

Applications

	Suggested dilution	Results
ELISA	1:20,000 - 1:60,000	
Western blot	0.2 µg/ml	Figure 1
Immunohistochemistry	1:100-1:500	
IF	1:100-1:500	

Target description

HDAC5 is a member of the class II mammalian histone deacetylase family, which is structurally related to yeast HDA1. Human HDAC5 is composed of 1122 amino acid residues. The deacetylase domain of HDAC5 is located at the C-terminal half of the molecule. The N-terminal non-deacetylase domain does not show any significant homology with any published sequence. Both domains are required for HDAC5-mediated repression of gene transcription. HDAC5 interacts with a growing number of transcriptional factors including MEF2A as well as other HDAC proteins. The interacting complexes bind to specific regions of chromatin and regulate gene transcription in these regions.

Results



Figure 1. HDAC5 antibody Western blot results

Western blot using the Diagenode HDAC5 antibody on mouse brain extracts (5 μ g per lane). The HDAC5 antibody was used at a dilution of 0.2 μ g/mL overnight at 4°C. Predicted/Observed size: 124 kDa.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing (Ougrée) - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: September 10, 2014