

## 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 µg per lane). The HDAC5 antibody was used at a dilution of 0.2 µg/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