

## Anti-HDAC7

Rabbit Polyclonal Antibody

**Catalog # H89-63R**

Lot # E208-8

### Cited Applications

Western blot (1:1000)

*Ideal working dilutions for each application should be empirically determined by the investigator.*

### Specificity

Recognizes the human HDAC7 protein

### Cross Reactivity

HDAC7 from other species may also be detectable

### Host

Rabbit

### Immunogen

Synthetic peptide corresponding to c-terminus of human HDAC7 conjugated to KLH.

### Formulation

PBS pH 7.4, 50% glycerol

### Stability

Store at 4°C (add 0.1% NaN<sub>3</sub>) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

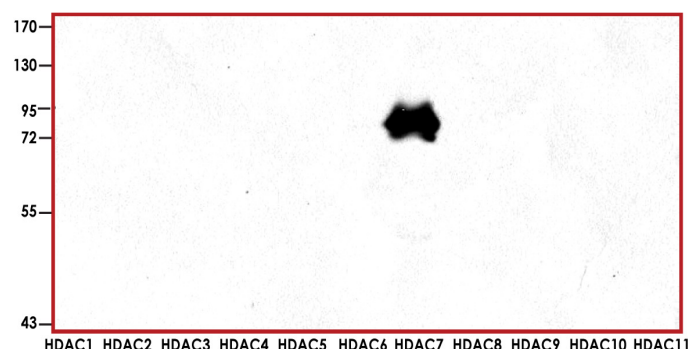
### Scientific Background

HDAC7 or Histone deacetylase 7 belongs to the histone deacetylase/acuc/apha family and is a component of the histone deacetylase complex. The protein encoded by HDAC7 gene has sequence homology to members of the histone deacetylase family whose protein promotes repression mediated via the transcriptional co-repressor SMRT (1). HDAC7 interacts with β-catenin keeping endothelial cells in a low proliferation stage. HDAC7 regulates NUR77 and apoptosis in developing thymocytes (2). HDAC7 is expressed in human osteoarthritis that contributes to cartilage degradation via promoting MMP-13 gene expression.

### References

1. Kao, H. Y. et.al: Isolation of a novel histone deacetylase reveals that class I and class II deacetylases promote SMRT-mediated repression. *Genes Dev.* 14: 55-66, 2000.
2. Dequiedt, F. et.al: HDAC7, a thymus-specific class II histone deacetylase, regulates Nur77 transcription and TCR-mediated apoptosis. *Immunity* 18: 687-698, 2003.

### Sample Data



Representative western blot with Anti-HDAC7 (1:1000) using 25 ng of human recombinant proteins HDAC1 to HDAC11.

## Anti-HDAC7

Rabbit Polyclonal Antibody

Catalog Number

H89-63R

Specific Lot Number

E208-8

Purification

Affinity chromatography

Concentration

0.25 µg/µL

Stability

1yr At -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**