

Anti-HDAC11

Rabbit Polyclonal Antibody

Catalog # H93-63R

Lot # E208-9

Cited Applications

Western blot (1:1000)

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

Specificity

Recognizes the human HDAC11 protein

Cross Reactivity

HDAC11 from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to c-terminus of human HDAC11 conjugated to KLH.

Formulation

PBS pH 7.4, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) 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.

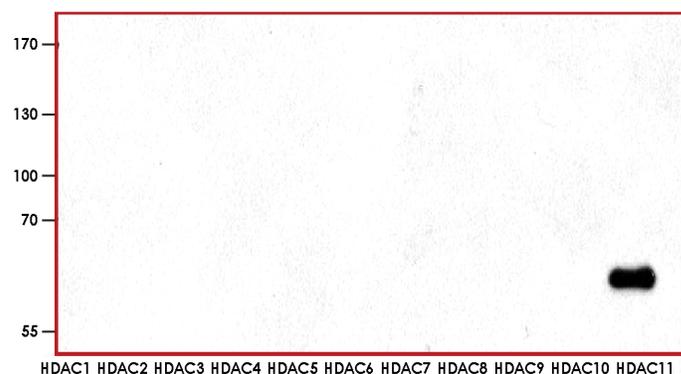
Scientific Background

HDAC11 belongs to the histone deacetylase/acuc/apha family and is a component of the histone deacetylase complex. HDAC11 belongs to class IV of the histone deacetylase family that localizes to the nucleus and is involved in regulating the expression of interleukin 10 (1). HDAC11 expression is the highest in brain, heart, skeletal muscle, kidney and testis. HDAC11 control the DNA expression by modifying the core histone octamers that package DNA into dense chromatin structures and repress gene expression. HDAC11 has HDAC activity and this activity is inhibitable by trapoxin, an HDAC inhibitor (2). In coimmunoprecipitation experiments, HDAC11 interacts with HDA6.

References

1. Villagra, A. et al: The histone deacetylase HDAC11 regulates the expression of interleukin 10 and immune tolerance. Nat Immunol. 2009 Jan;10(1):92-100.
2. Gao, L. et.al: Cloning and functional characterization of HDAC11, a novel member of the human histone deacetylase family. J. Biol. Chem. 277: 25748-25755, 2002.

Sample Data



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

Anti-HDAC11

Rabbit Polyclonal Antibody

Catalog Number

H93-63R

Specific Lot Number

E208-9

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
www.signalchem.com

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