

Anti-HDAC5

Rabbit Polyclonal Antibody

Catalog # H87-63R

Lot # E208-6

Cited Applications

Western blot (1:1000)

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

Specificity

Recognizes the human HDAC5 protein

Cross Reactivity

Only a weak reaction with human HDAC4 in Western Blot.

HDAC5 from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to c-terminus of human HDAC5 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.

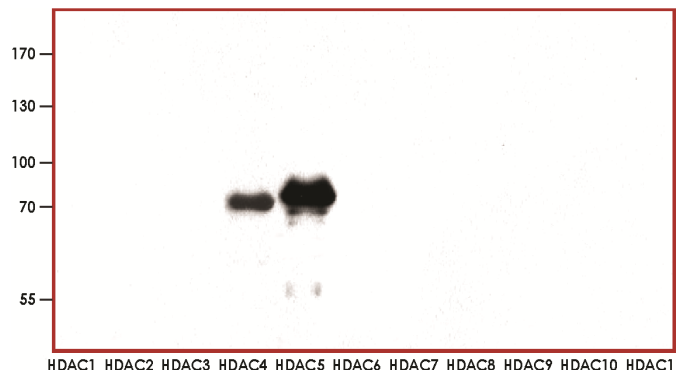
Scientific Background

HDAC5 or Histone deacetylase 5 belongs to the class II histone deacetylase/acuc/apha family that possesses histone deacetylase activity and represses transcription when tethered to a promoter (1). HDAC 5 plays a critical role in transcriptional regulation, cell cycle progression, and developmental events and also acts as a potential therapeutic target for the prevention of atherosclerosis. HDAC5 can co-immunoprecipitates with HDAC3 family members forming multicomplex proteins. HDAC5 can also interact with myocyte enhancer factor-2 (MEF2) proteins, resulting in repression of MEF2-dependent genes. HDAC5 gene is thought to be associated with colon cancer (2).

References

1. Grozinger, C. M. et.al : Three proteins define a class of human histone deacetylases related to yeast Hda1p. Proc. Nat. Acad. Sci. 96: 4868-4873, 1999.
2. Scanlan, M. J. et.al: Characterization of human colon cancer antigens recognized by autologous antibodies. Int. J. Cancer 76: 652-658, 1998.

Sample Data



Representative western blot with Anti-HDAC5 (1:1000) using 25 ng of human recombinant proteins HDAC1 to HDAC11.

Anti-HDAC5

Rabbit Polyclonal Antibody

Catalog Number

H87-63R

Specific Lot Number

E208-6

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.