

Anti-HDAC6

Rabbit Polyclonal Antibody

Catalog # H88-63R

Lot # E208-7

Cited Applications

Western blot (1:1000)

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

Specificity

Recognizes the human HDAC6 protein

Cross Reactivity

Weak reaction with human HDAC11 in Western Blot.

HDAC6 from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to c-terminus of human HDAC6 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.

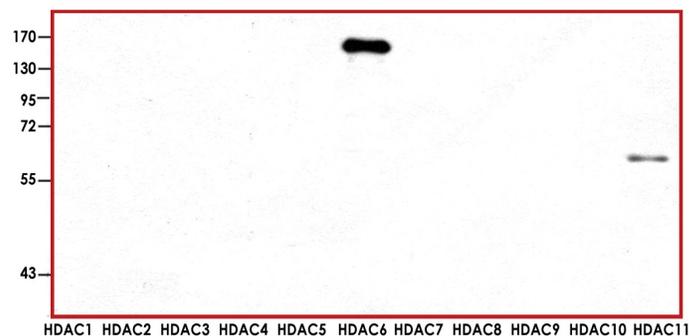
Scientific Background

HDAC6 or Histone deacetylase 6 belongs to the histone deacetylase/acuc/apha family and is a component of the histone deacetylase complex. Histones play a critical role in transcriptional regulation, cell cycle progression, and developmental events. The protein encoded by this gene belongs to class II of the histone deacetylase family that regulates important biologic processes beyond histone metabolism and gene transcription (1). HDAC6 is a central component of the stress response that regulates SG formation and potentially contributes to control of RNA metabolism and translation (2).

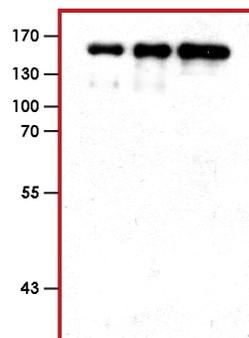
References

1. Hubbert, C. et.al: HDAC6 is a microtubule-associated deacetylase. Nature 417: 455-458, 2002.
2. Kwon, S. et.al: The deacetylase HDAC6 is a novel critical component of stress granules involved in the stress response. Genes Dev. 21: 3381-3394, 2007.

Sample Data



Representative western blot with Anti-HDAC6 (1:1000) using 50ng of human recombinant proteins HDAC1 to HDAC11.



Representative western blot with Anti-HDAC6 (1:1000) using 12ng, 25ng and 50ng of recombinant GST-HDAC6 protein.

Anti-HDAC6

Rabbit Polyclonal Antibody

Catalog Number

H88-63R

Specific Lot Number

E208-7

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.