

Anti-phospho-HDAC2 (Ser394)

Rabbit Polyclonal Antibody

Catalog # H84-65R

Lot # Z2014-37

Cited Applications

IHC, IF, WB

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

Specificity

Recognizes the HDAC2 protein phosphorylated at serine 394

Cross Reactivity

Human, Mouse, and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Phosphopeptide derived from human HDAC2 around the phosphorylation site of serine 394 (E-D-SP-G-D).

Formulation

Phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

Stability

1yr at -20°C from date of shipment

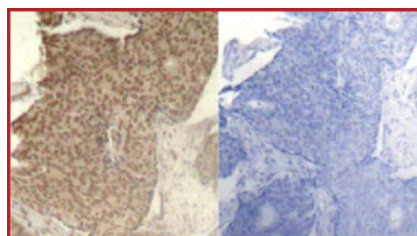
Scientific Background

HDAC2 or Histone deacetylase 2 belongs to the histone deacetylase family that acts via the formation of large multiprotein complexes, and is responsible for the deacetylation of lysine residues at the N-terminal regions of core histones (H2A, H2B, H3 and H4). HDAC2 forms transcriptional repressor complexes by associating with many different proteins and plays an important role in transcriptional regulation, cell cycle progression and developmental events. HDAC2 functions in modulating synaptic plasticity and long-lasting changes of neural circuits, which in turn negatively regulates learning and memory (1). HDAC1 and HDAC2 are functionally redundant in cardiac growth and development and they maintain cardiomyocyte identity and function (2).

References

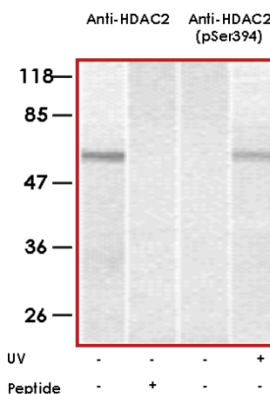
- Guan, J.S. et.al: HDAC2 negatively regulates memory formation and synaptic plasticity. *Nature* 459: 55-60, 2009.
- Montgomery, R. L. et.al: Histone deacetylases 1 and 2 redundantly regulate cardiac morphogenesis, growth, and contractility. *Genes Dev.* 21: 1790-1802, 2007.

Sample Data



P-Peptide - +

Immunohistochemical analysis of paraffin-embedded human breast carcinoma tissue using Anti-phospho-HDAC2 (Ser394).



Western blot analysis of extracts from HT-29 cells untreated or treated with UV (20 mins), using HDAC2 antibody and Anti-phospho-HDAC2 (Ser394) (Lane 3 and 4)

Anti-phospho-HDAC2 (Ser394)

Rabbit Polyclonal Antibody

Catalog #

H84-65R

Lot #

Z2014-37

Purification

Affinity chromatography

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.