

HDAC11 antibody (pAb)

Catalog No: 39208**RRID:** AB_2614938**Application(s):** WB**Reactivity:** Human, Mouse, Rat**Volume:** 200 µl**Purification:** Affinity Purified**Host:** Rabbit**Isotype:** IgG**Concentration:** 0.5 µg/µl**Molecular Weight:** 36 kDa

Background: HDAC11 (Histone Deacetylase 11) is a member of the class IV mammalian histone deacetylases (HDACs) involved in regulating chromatin structure during transcription. These enzymes catalyze the removal of acetyl groups from lysine residues of histones and other cellular proteins. Lysine N-ε-acetylation is a dynamic, reversible and tightly regulated protein and histone modification that plays a major role in regulation of gene expression in various cellular functions. It consists of the transfer of an acetyl moiety from an acetyl coenzyme A to the ε-amino group of a lysine residue.

In vivo, acetylation is controlled by the antagonistic activities of histone acetyltransferases (HATs) and histone deacetylases (HDACs). The HDACs are grouped into four classes, on the basis of similarity to yeast counterparts: HDAC class I (HDAC1, HDAC2, HDAC3 and HDAC8), class II (HDAC4, HDAC5, HDAC6, HDAC7, 9 and 10), class III (SIRT1-7) and class IV (HDAC11). In humans, expression of HDAC11 is limited to only a few organs, including brain, heart, muscle and kidney, suggesting an important role of HDAC11 in those tissues.

Immunogen: This HDAC11 antibody was raised against a peptide corresponding to amino acid residues 238-251 of human HDAC11.

Buffer: PBS containing 0.2% gelatin and 0.05% sodium azide. Sodium azide is highly toxic.

Application Notes:

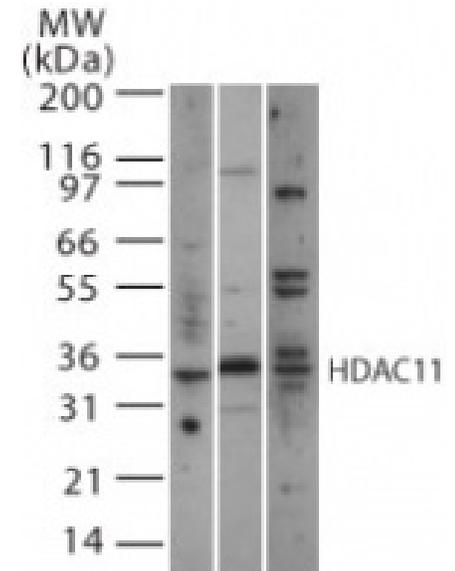
Applications Validated by Active Motif:

WB: 1 - 3 µg/ml dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Store at 4°C for short term. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

**HDAC11 pAb tested by Western blot.**

Detection of HDAC11 by Western blot. The analysis was performed using HDAC11 pAb at a dilution of 2 µg/ml and brain whole-cell extracts from human (lane 1), mouse (lane 2) or rat origin.