

Histone macroH2A2 antibody (pAb)

Catalog Nos: 39873, 39874

RRID: AB_2793376

Isotype: IgG

Application(s): ICC, IF, IHC, WB

Reactivity: Human

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 40 kDa

Background: Histone macroH2A2 is a histone H2A variant and one of the three members of the macroH2A family. The three macroH2A proteins contain the C-terminal 'macro' domain in addition to the H2A homologous region. This macro domain may bind poly ADP-ribose and other NAD⁺ metabolites. The macroH2A1.1 and macroH2A1.2 proteins are produced from alternatively spliced transcripts from the macroH2A1 gene while a second distinct macroH2A gene encodes macroH2A2. In mammals, macroH2A is preferentially localized to the inactive X chromosome and in general is proposed to play a role in repressing gene expression.

Immunogen: This macroH2A2 antibody was raised against a peptide corresponding to amino acids 162-372 of human macroH2A2.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

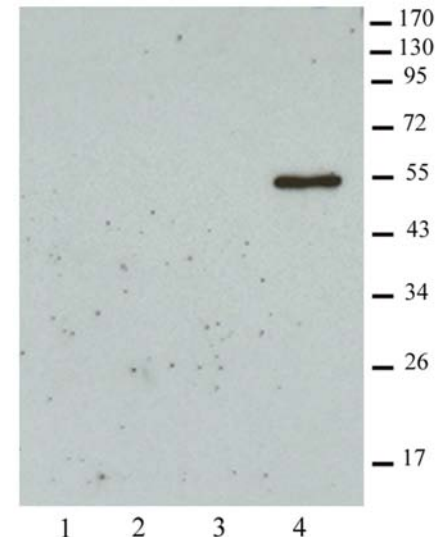
Application Notes:

Applications Validated by Active Motif:

WB: 1:500 - 1:1,000 dilution

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. 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.



Histone macroH2A2 antibody (pAb) tested by Western blot.

U2OS cells mock transfected (lane 1) or transfected with GST-Tagged macroH2A1.1 (lane 2), GST-macroH2A1.2 (lane 3) or GST-macroH2A2 (lane 4) were probed with Histone macroH2A2 antibody at a dilution of 1:1,000.