

## Histone macroH2A1.2 antibody (mAb)

**Catalog Nos:** 61427, 61428

**RRID:** AB\_2793631

**Clone:** 14G7-H9

**Isotype:** IgG2b

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

**Reactivity:** Human

**Volumes:** 100  $\mu$ l, 10  $\mu$ l

**Purification:** Ascites

**Host:** Mouse

**Molecular Weight:** 40 kDa

**Background:** Histone macroH2A1.2 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<sup>+</sup> 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. In mice, macroH2A1 is required for the silencing of endogenous retroviruses.

**Immunogen:** Recombinant protein corresponding to amino acids 198-228 of human Histone macroH2A1.2.

**Buffer:** Ascites containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

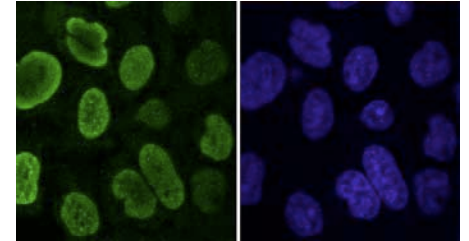
Applications Validated by Active Motif:

WB: 1:300 - 1:1,000 dilution

ICC/IF: 1:50 - 1:200 dilution

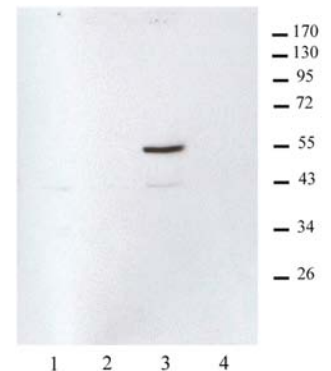
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This 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 macroH2A1.2 antibody (mAb) tested by immunofluorescence.

Staining of macroH2A1.2 transfected HeLa cells with Histone macroH2A1.2 antibody (1:50 dilution, left panel) and DAPI (right panel).



### Histone macroH2A1.2 antibody (mAb) tested by Western blot.

U2OS cells mock transfected (lane 1) or transfected with tagged macroH2A1.1 (lane 2), macroH2A1.2 (lane 3) or macroH2A2 (lane 4) were probed with Histone macroH2A1.2 antibody at a dilution of 1:300. Band at 40kD is the endogenous macroH2A1.2.