

Histone H2A.Z.2.2 antibody (mAb)

Catalog Nos: 61737, 61738

RRID: AB_2793750

Clone: 1H11-11

Isotype: IgG1

Application(s): WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 12 kDa

Background: **Histone H2A.Z** (H2A Histone Family, Member V) is a histone H2A variant, a protein similar to canonical H2A but with different molecular identity and unique functions. H2A.Z is highly conserved during evolution. It plays an important role in basic cellular mechanisms such as gene activation, chromosome segregation, heterochromatic silencing and progression through the cell cycle. H2A.Z is acetylated at multiple lysine residues in its amino terminus, which may serve to allow H2A.Z to function as an insulator of chromatin domains. In the case of **H2A.Z.2.2**, this variant has been shown to be responsible for destabilization of the nucleosome through the chaperone complexes TIP60 and SRCAP.

Immunogen: This antibody was raised against a peptide within the C-terminal region of human Histone H2A.Z.2.2.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

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.