

HIRA antibody (mAb)

Catalog Nos: 39555, 39556

RRID: AB_2793255

Clone: WC15

Isotype: IgG

Application(s): IP, WB

Reactivity: Human, Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 120 kDa

Background: HIRA (DGCR1, TUP1) is a histone chaperone that preferentially places the variant Histone H3.3 in nucleosomes, replacing canonical Histone H3 during the cell cycle. HIRA is required for the periodic repression of histone gene transcription during the cell cycle and when complexed with ASF1α is involved in the formation of the senescence-associated heterochromatin foci.

Immunogen: This HIRA antibody was raised against a GST-fusion protein containing residues 421-729 of human HIRA.

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:

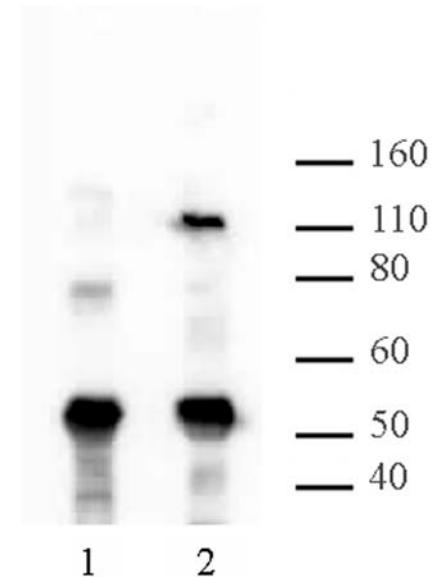
Applications Validated by Active Motif:

IP: 5 µg per IP

WB: 2 - 5 µg/ml 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.



HIRA mAb (Clone WC15) tested by Western blot.

5 µg of HIRA mAb (Clone WC15) was used to immunoprecipitate HIRA from 500 µg of colcemid-treated U2OS nuclear extract (lane 2). 5 µg of control mouse IgG was also used (lane 1). The immunoprecipitated protein was detected by Western blotting using HIRA mAb (Clone WC119.2H11) at a dilution of 5 µg/ml.