

## DHX9 antibody (mAb)

**Catalog Nos:** 61705, 61706

**RRID:** AB\_2793743

**Clone:** 8E3

**Isotype:** IgG1

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

**Reactivity:** Human, Mouse

**Quantities:** 100 µg, 10 µg

**Purification:** Protein A Chromatography

**Host:** Rat

**Concentration:** 1 µg/µl

**Molecular Weight:** 150 kDa

**Background:** DHX9 (DEAH (Asp-Glu-Ala-His) Box Helicase 9) unwinds double-stranded DNA and RNA in a 3' to 5' direction. Alteration of secondary structure may subsequently influence interactions with proteins or other nucleic acids. Functions as a transcriptional activator. Component of the CRD-mediated complex that promotes MYC mRNA stability. Involved with LARP6 in the stabilization of type I collagen mRNAs for CO1A1 and CO1A2. Positively regulates HIV-1 LTR-directed gene expression.

**Immunogen:** This antibody was raised against a peptide within the N-terminal region of mouse DHX9.

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

### Application Notes:

Applications Validated by Active Motif:

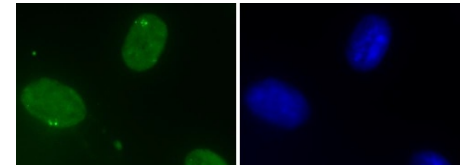
ICC/IF: 2 - 10 µg/ml dilution

WB: 0.5 - 2 µg/ml dilution

The addition of 0.1% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

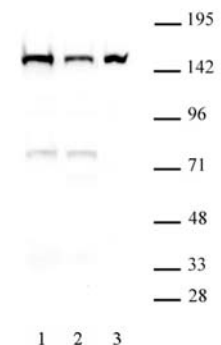
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This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



### DHX9 antibody (mAb) (Clone 8E3) tested by Immunofluorescence.

Left: Formaldehyde fixed L929 cells stained with DHX9 antibody (mAb). Right: Hoechst.



### DHX9 antibody (mAb) (Clone 8E3) tested by Western blot.

DHX9 antibody detection by Western blot. The analysis was performed using 30 µg of either NIH-3T3 nuclear extract (lane 1), NIH-3T3 cytoplasmic extract (lane 2), or HeLa nuclear extract (lane 3) and DHX9 antibody (mAb) at a 2 µg/ml dilution.