

MEIS 1/2/3 antibody (mAb)

Catalog Nos: 39795, 39096, 39796

RRID: AB_2750570

Clone: 9.2.7

Isotype: IgG1

Application(s): ICC, IF

Reactivity: Human, Mouse

Quantities: 100 µg, 50 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 54-66 kDa

Background: The MEIS family of proteins (Myeloid Ecotropic viral Integration Site homologues) encode proteins that function as cofactors for the Homeobox (HOX) proteins. Homeobox genes play a crucial role in normal development. In addition, several homeoproteins are involved in neoplasia. MEIS and a functionally related protein Pbx belong to the TALE (Three Amino acid Loop Extension) homeodomain superfamily of proteins. The role of MEIS proteins in the regulation of transcription is as yet unclear.

Immunogen: This MEIS antibody was raised against a recombinant protein corresponding to amino acids 50-390 mouse MEIS1.

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:

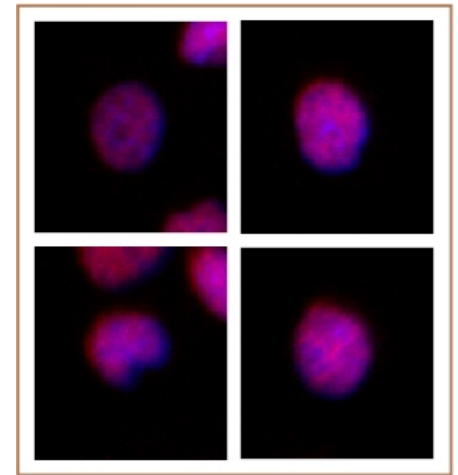
Applications Validated by Active Motif:

ICC/IF: 0.5 µg/ml dilution

This antibody is also available as an AbFlex® engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91259.

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.



MEIS 1/2/3 antibody (mAb) tested by immunofluorescence.

Formaldehyde fixed HeLa cells stained with MEIS 1/2/3 antibody at a 0.5 µg/ml dilution.