

## MAZ antibody (mAb)

**Catalog Nos:** 39935, 39936

**RRID:** AB\_2793398

**Clone:** 133

**Isotype:** IgG

**Application(s):** ChIP, EMSA, WB

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 55 kDa

**Background:** MAZ (Myc-associated zinc finger protein) functions as a transcription factor roles in both transcription initiation and termination. MAZ binds to two sites within the c-Myc promoter, ME1a1 and ME1a2. It also binds to multiple G/C-rich sites within the promoter of the Sp1 family of transcription factors. MAZ also participates in skeletal muscle development through the activation of several skeletal muscle-specific genes. MAZ has also been reported to drive tumor-specific expression of PPAR-gamma in breast cancer cells.

**Immunogen:** This MAZ antibody was raised against recombinant human MAZ protein.

**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.

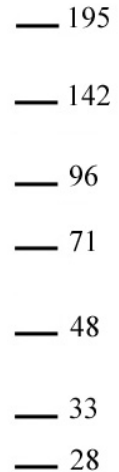
### Application Notes:

Applications Validated by Active Motif:  
WB

This antibody is also available as an AbFlex<sup>®</sup> engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91205.

**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.



**MAZ antibody (mAb) tested by Western Blot.**  
20 µg MCF7 cell nuclear extract was run on SDS-PAGE, and probed with 2 µg/ml MAZ antibody.  
MW: 55 kDa