

## Mi-2 beta antibody (mAb)

**Catalog Nos:** 39695, 39696

**RRID:** AB\_2793309

**Clone:** 3F2/4

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human, Mouse

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 240 kDa

**Background:** Mi-2β (Mi-2beta, CHD4, chromodomain helicase DNA-binding protein 4) represents the major Mi-2 isoform present in the complex Mi-2/NuRD. This complex is a multi-subunit protein complex containing both histone deacetylase and nucleosome-dependent ATPase subunits. Current models predict that this complex functions primarily in transcriptional repression. In addition to the ATPase domain, Mi-2 beta contains a chromodomain and two PHD domains. Mi-2 beta interacts with BRG-1, a component of the SWI/SNF complex, and the RET finger protein (RFP).

**Immunogen:** This Mi-2 beta antibody was raised against a synthetic peptide derived from human Mi-2B.

**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:

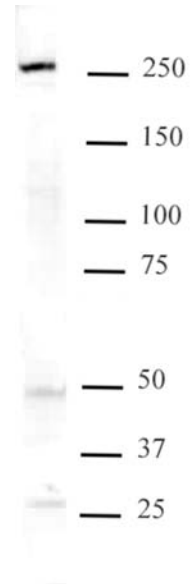
Applications Validated by Active Motif:

WB: 1 - 2 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. 91275.

**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.



### Mi-2 beta pAb tested by Western blot.

Nuclear extract of Jurkat cells stained with Mi-2 beta mAb at a dilution of 2 µg/ml.