

MXD1 antibody (mAb)

Catalog Nos: 61739, 61740

RRID: AB_2793751

Clone: 5F4

Isotype: IgG2b

Application(s): ChIP, IP, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 30 kDa

Background: **MXD1** (MAX Dimerization Protein 1) is a transcriptional repressor. MAD binds with MAX to form a sequence-specific DNA-binding protein complex which recognizes the core sequence 5-CAC[GA]TG-3. MAD thus antagonizes MYC transcriptional activity by competing for MAX.

Immunogen: This antibody was raised against a recombinant protein corresponding amino acids 6-221 of human MXD1.

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

Application Notes:

Published Applications:

ChIP

IP

WB

See reference for more information. 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.