

MRG15 antibody (pAb)

Catalog Nos: 39361, 39362

RRID: AB_2615022

Isotype: IgG

Application(s): WB

Reactivity: Human, Wide Range Predicted

Volumes: 200 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 39 kDa

Background: MRG15 (MORF4-Related Gene on chromosome 15) is the homolog of the yeast Eaf3 protein and belongs to the MRG family of genes, which are involved in various cellular processes as cellular aging, positive and negative transcriptional regulation and DNA damage repair. MRG15 can be associated with both histone deacetylase and histone acetyltransferase complexes (such as the NuA4 histone acetyltransferase complex, which is involved in transcriptional activation of select genes principally by acetylation of nucleosomal Histone H2A and Histone H4). MRG15, thus, potentially plays a role in both transcriptional silencing and activation. In addition to the more conserved MRG motifs, MRG15 also contains a chromodomain near the N-terminus.

Immunogen: This MRG15 antibody was raised against a peptide corresponding to the C-terminus of human MRG15.

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.

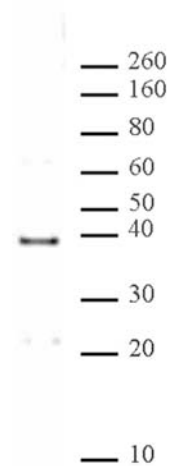
Application Notes:

Applications Validated by Active Motif:

WB: 1:250 - 1:1,000 dilution

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.



MRG15 pAb tested by Western blot.

Nuclear extract of HeLa cells (20 µg) was probed with MRG15 pAb (1:500 dilution).