

RYPB antibody (pAb)

Catalog Nos: 61461, 61462

RRID: AB_2793647

Isotype: IgG

Application(s): WB

Reactivity: Human, Mouse

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 35 kDa

Background: RYPB (RING1 and YY1 Binding Protein) is a member of a non-canonical PRC1 repressive complex composed of at least RING1B and ubiquitinated H2A. Also inhibits ubiquitination and subsequent degradation of TP53, and thereby plays a role in regulating transcription of TP53 target genes. May be implicated in the regulation of transcription as a repressor of the transcriptional activity of E4TF1.

Immunogen: This antibody was raised against a peptide within the C-terminal region of human RYPB.

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

Application Notes:

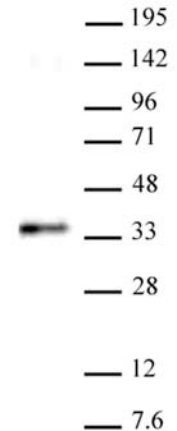
Applications Validated by Active Motif:

WB*: 1:500 - 1:2,000 dilution

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

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.



RYPB antibody (pAb) tested by Western blot.
Nuclear extract of mouse ES cells (30 μ g) probed with RYPB antibody (pAb) at a dilution of 1:500.