

PHB / Prohibitin antibody (pAb)

Catalog Nos: 61217, 61218

RRID: AB_2793555

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 30 kDa

Background: PHB / Prohibitin is a key cell cycle regulatory protein, inhibiting cell division by inactivating E2F family proteins, and has potent anti-proliferative activity. PHB / Prohibitin is also involved in repression of transcription through androgen receptor.

Immunogen: This PHB / prohibitin antibody was raised against a peptide within the N-terminal region of human PHB.

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

Application Notes:

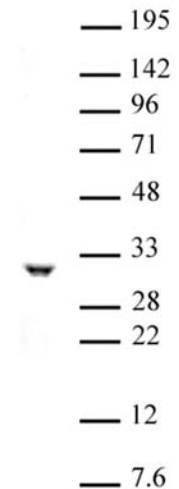
Applications Validated by Active Motif:

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

For optimal results, primary antibody incubations should be performed at 4°C. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. 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.



PHB / Prohibitin antibody (pAb) tested by Western blot.

Whole-cell extract of MCF-7 cells (25 µg) probed with PHB antibody (pAb) at a dilution of 1:500.