

Rad18 antibody (mAb)

Catalog No: 40979

RRID: AB_2793469

Clone: 79B1048

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Mouse

Isotype: IgG1

Concentration: 0.5 µg/µl

Molecular Weight: 60 kDa

Background: Rad18 (RNF73, hHR18) is highly similar to *S. cerevisiae* DNA damage repair protein Rad18. Yeast Rad18 functions through its interaction with Rad6, which is an ubiquitin-conjugating enzyme required for post-replication repair of damaged DNA. Rad18 is similar to its yeast counterpart, as it is able to interact with the human homolog of yeast Rad6 protein through a conserved ring-finger motif.

Immunogen: This Rad18 antibody was raised against a synthetic peptide corresponding to amino acid residues 402-414 of human Rad18.

Buffer: PBS containing 0.02% sodium azide. Sodium azide is highly toxic.

Application Notes:

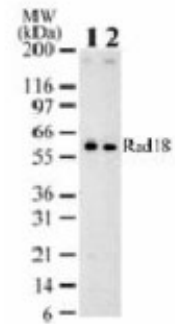
Applications Validated by Active Motif:

WB: 1 - 2 µg/ml dilution

For optimal results, primary antibody incubations should be performed at room temperature. 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. Store at 4°C for short term. 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.

**Rad18 mAb tested by Western blot.**

Rad18 detection by Western blot. hRad18 mAb was diluted at 2 µg/ml (lane 1) and 1 µg/ml (lane 2) and tested against 10 µg of HL-60 nuclear extract by Western blot. A single band at 60 kDa is detected.