

## Rad17 phospho Ser647 antibody (pAb)

**Catalog No:** 39222

**RRID:** AB\_2793192

**Application(s):** WB

**Reactivity:** Human, Rat

**Quantity:** 100 µg

**Purification:** Affinity Purified

**Host:** Rabbit

**Isotype:** IgG

**Concentration:** 0.5 µg/µl

**Molecular Weight:** 77 kDa

**Background:** Rad17 Ser647 (Rad24, CCYC, R24L) is essential for sustained cell growth, maintenance of chromosomal stability, and ATR-dependent checkpoint activation upon DNA damage. Rad17 has a weak ATPase activity required for binding to chromatin. Rad17 participates in the recruitment of the Rad1-RAD9-HUS1 complex onto chromatin, and in CHEK1 activation. Rad17 may also serve as a sensor of DNA replication progression, and may be involved in homologous recombination.

**Immunogen:** This Rad17 phospho Ser647 antibody was raised against a synthetic peptide corresponding to mouse Rad17 phosphorylated at serine 647 .

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

### Application Notes:

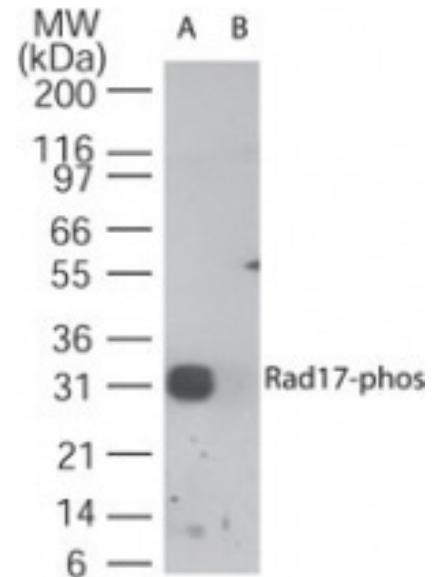
Applications Validated by Active Motif:

WB: 0.1 - 1 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.



### Rad17 phospho Ser647 pAb tested by Western blot.

Rad17 detection by Western blot. The analysis of phospho-RAD17 was performed using the Rad17 phospho Ser647 pAb and partial recombinant Rad17 protein containing phospho Ser647 (lane A) and the same sequence without a phospho serine (lane B).