

## Rad21 antibody (mAb)

**Catalog Nos:** 39383, 39384

**RRID:** AB\_2688003

**Clone:** 52A311

**Isotype:** IgG

**Application(s):** ICC, IF, WB

**Reactivity:** Human, Mouse

**Quantities:** 200 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 100 kDa

**Background:** Rad21 Clone 52A311 (HR21, MCD1, NXP1, hHR21) is highly similar to the gene product of *S. pombe* Rad21, a gene involved in the repair of DNA double-strand breaks, as well as in chromatid cohesion during mitosis. Rad21 is a nuclear phospho-protein, which becomes hyperphosphorylated in cell cycle M phase. The highly regulated association of Rad21 with mitotic chromatin specifically at the centromere region suggests its role in sister chromatid cohesion in mitotic cells.

**Immunogen:** This Rad21 antibody was raised against recombinant protein corresponding to the N-terminal half of mouse Rad21.

**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:

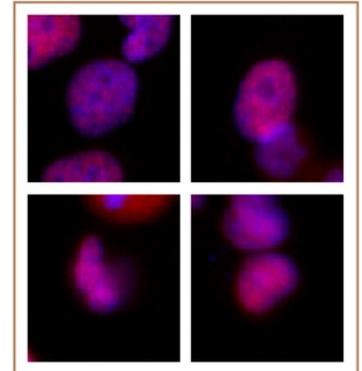
ICC/IF: 0.5 µg/ml dilution

WB: 0.5 - 2 µg/ml dilution

This antibody is also available as an AbFlex<sup>®</sup> engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91245.

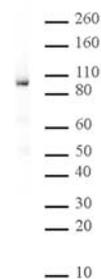
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This 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.



### Rad21 antibody (mAb) tested by immunofluorescence.

Formaldehyde fixed HeLa cells stained with Rad21 antibody at a 0.5 µg/ml dilution.



### Rad21 antibody (mAb) tested by Western blot.

HeLa nuclear extract (30 µg per lane) was probed with Rad21 mAb (Clone 52A311) (1 µg/ml).