

Rad51 antibody (pAb)

Catalog Nos: 39194, 39195

RRID: AB_2614951

Isotype: IgG

Application(s): ICC, IF

Reactivity: Human

Quantities: 200 µg, 10 µg

Purification: Protein A Chromatography

Host: Rabbit

Concentration: 1 µg/µl

Molecular Weight: 37 kDa

Background: Rad51 (BRCC5, RecA) is known to be involved in the homologous recombination and repair of DNA. Rad51 can interact with the ssDNA-binding protein RPA and Rad52, and it is thought to play roles in homologous pairing and strand transfer of DNA. Rad51 is also found to interact with BRCA1 and BRCA2, which may be important for the cellular response to DNA damage. BRCA2 is shown to regulate both the intracellular localization and DNA-binding ability of this protein. Loss of these controls following BRCA2 inactivation may be a key event leading to genomic instability and tumorigenesis.

Immunogen: This Rad51 antibody was raised against full-length human Rad51 expressed in *E. coli*.

Buffer: PBS pH 7.5 containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

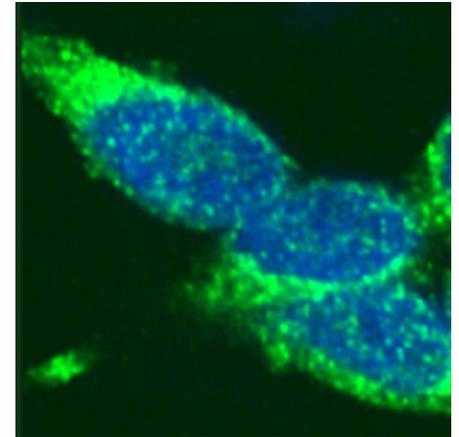
Application Notes:

Applications Validated by Active Motif:

ICC/IF: 1 - 5 µg/ml dilution

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.



Rad51 pAb tested by immunofluorescence. HEK293 cells stained with Rad51 polyclonal antibody (green) and DAPI (blue).