

Rad52 antibody (pAb)

Catalog Nos: 39196, 39197

RRID: AB_2793179

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: 46 kDa

Background: Rad52 is important for DNA double-strand break repair and homologous recombination. Rad52 is shown to bind single-stranded DNA ends, and mediate the DNA-DNA interaction necessary for the annealing of complementary DNA strands. Rad52 is also found to interact with DNA recombination protein Rad51, which suggested its role in RAD51-related DNA recombination and repair.

Immunogen: This Rad52 antibody was raised against full-length human Rad52 expressed in *E. coli*.

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

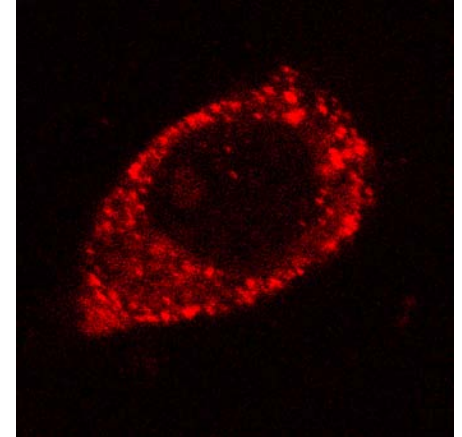
Application Notes:

Applications Validated by Active Motif:

ICC/IF: 2 - 10 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.



Rad52 pAb tested by immunofluorescence.
Staining of HEK293 cells with Rad52 pAb at a 2 µg/ml dilution.