



Anti-XRCC2

CATALOG NO.: 53956

BACKGROUND:

XRCC2 (X-Ray Cross Complementing 2) is a 34 kDa protein that is essential for repairing the double strand DNA breaks by homologous recombination between sister chromatids. It is homologous to Rad51. Five human homologues of RAD 51; XRCC2, XRCC3, RAD51B, RAD51C, RAD51D, interact among themselves to form one big complex or several smaller complexes.

SOURCE & REACTIVITY:

Rabbit anti-XRCC2 polyclonal antibody was raised against a synthetic peptide derived from the N-terminus of human XRCC2. Species reactivity includes human, while others are not tested.

APPLICATION:

The following concentration ranges are recommended starting points for this product.

WB

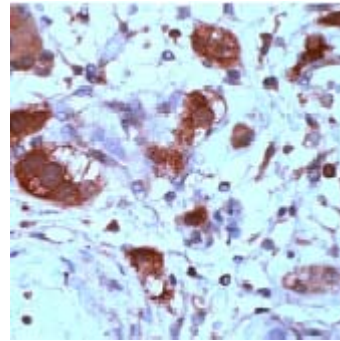
IHC (Formalin/paraffin)

Working dilution: 1:200 for 10 min at room temperature (RT)

Staining of formalin-fixed tissues requires boiling tissue sections in 10mM citrate buffer, pH 6.0 for 10 min followed by cooling at RT for 20 min.

Positive Control: Testis

Staining Pattern: Cytoplasmic, nuclear



Human testis
stained with
anti-XRCC2
(Cat. #53956)

This product is for in vitro research purposes only.

STORAGE:

0.25 ml of affinity purified antibody at 0.2 mg/ml in 1X PBS (pH 7.4) with 0.2% BSA and 0.05% sodium azide. Store at 2-8 °C for up to 1 year. Avoid repeated freeze thaw cycles.