

WRN antibody (mAb)

Catalog Nos: 61169, 61170

RRID: AB_2793536

Clone: 195C

Isotype: IgG1

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

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 162 kDa

Background: WRN protein is a RECQ helicase that contains 3' to 5' helicase activity and DNA-dependent ATPase activity. WRN takes part in many processes, including DNA replication, recombination, repair, and telomere maintenance. Mutations in the gene encoding WRN cause an autosomal recessive segmental progeria called **Werner syndrome** (WS). WS patients prematurely develop atrophic skin, thin gray hair, osteoporosis, type II diabetes, cataracts, arteriosclerosis and cancer.

Immunogen: This WRN antibody was raised against a recombinant protein corresponding to amino acids 1072 - 1432 of human WRN.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

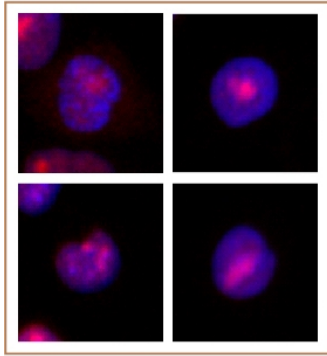
ICC/IF: 0.5 µg/ml dilution

IHC(P): User optimization required

WB: 0.5 - 2 µg/ml dilution

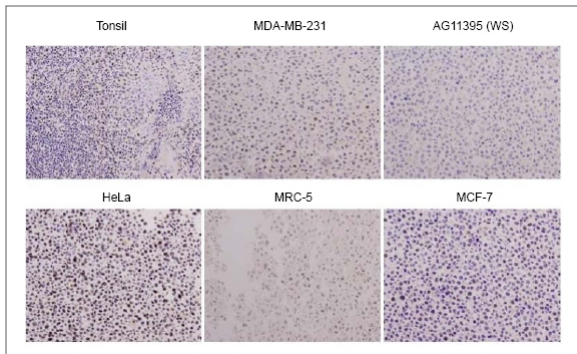
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



WRN antibody (mAb) tested by Immunofluorescence.

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



WRN antibody (mAb) tested by Immunohistochemistry.

Paraffin embedded tissue samples: human tonsil, MDA-MB-231 (breast adenocarcinoma cell line), AG11395 WS (fibroblast of a patient with Werner Syndrome immortalized with SV40), HeLa cells, MRC-5 (lung fibroblast cell line), and MCF-7 (breast carcinoma cell line).



WRN antibody (mAb) tested by Western blot.

Detection of WRN by Western blot. Lane 1: HeLa cell extract. Lane 2: WS cell extract (fibroblast from a patient with Werner Syndrome).