

## ATM phospho Ser1981 antibody (mAb)

**Catalog Nos:** 39529, 39530

**RRID:** AB\_2793244

**Clone:** 10H11.E.12

**Isotype:** IgG1, k

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

**Reactivity:** Human, Mouse

**Volumes:** 200  $\mu$ l, 10  $\mu$ l

**Purification:** Ascites

**Host:** Mouse

**Molecular Weight:** 370 kDa

**Background:** ATM – Ataxia Telangiectasia Mutated is a PI3-family protein kinase and a critical cell cycle checkpoint protein. ATM and a related protein, ATR, play crucial roles in the maintenance of genome integrity and the response to DNA damage. Upon activation, ATM phosphorylates itself at serine 1981 and then phosphorylates a number of downstream proteins (*e.g.* BRCA1, CHK2, H2AX, p53, NBS1 and SMC1) resulting in cell cycle arrest and the initiation of DNA damage repair. Loss of function of ATM is causal to a variety of syndromes involving increased incidence of several cancers.

**Immunogen:** This ATM phospho Ser1981 antibody was raised against a peptide containing phospho-serine 1981 of human ATM.

**Buffer:** Ascites containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

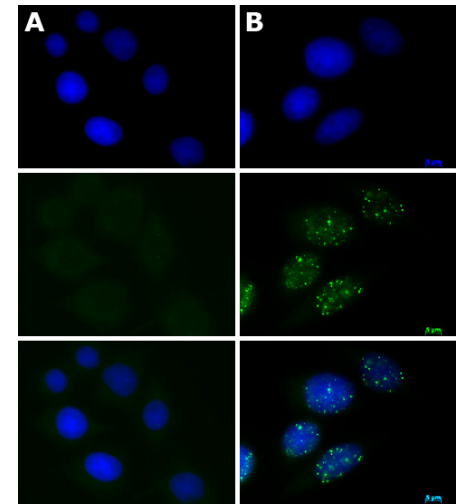
Applications Validated by Active Motif:

ICC/IF: 1:500 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. 91207.

**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.



### ATM phospho Ser1981 antibody (Clone 10H11.E.12) tested by immunofluorescence.

HeLa cells stained with ATM phospho Ser1981 antibody (Clone 10H11.E.12) at a 1:500 dilution using using MAX Stain<sup>™</sup> Immunofluorescence Tools.

Panel A: Normal HeLa cells.

Panel B: HeLa cells treated with 3 Gy ionizing radiation and collected 30 minutes post-exposure.

Top images: Cells stained with DAPI.

Middle images: Same cells stained with ATM phospho Ser1981 mAb (Clone 10H11).