

## MOF / MYST1 antibody (pAb)

**Catalog Nos:** 61245, 61246

**RRID:** AB\_2793566

**Isotype:** IgG

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

**Reactivity:** Human

**Volumes:** 100  $\mu$ l, 10  $\mu$ l

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 65 kDa

**Background:** **MOF / MYST1** (ortholog of *Drosophila* males absent on the first (MOF), MYST histone acetyltransferase 1, Histone acetyltransferase KAT8) is a histone acetyltransferase which may be involved in transcriptional activation. Member of the MYST family of histone acetyltransferases. Contains a chromodomain and C2HC-type zinc finger motif. Component of a multisubunit histone acetyltransferase complex (MSL) at least composed of the MOF/KAT8, MSL1/hampin, MSL2L1 and MSL3L1. Interacts with ATM through the chromodomain and may influence the function of ATM.

**Immunogen:** This MOF / MYST1 antibody was raised against a peptide within the N-terminal region of human MOF / MYST1 (MYST histone acetyltransferase 1).

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

### Application Notes:

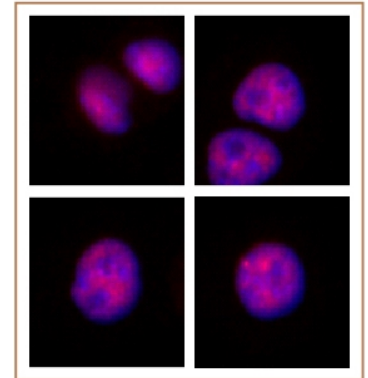
Applications Validated by Active Motif:

ICC/IF: 1:2,000 dilution

WB: 1:500 - 1:1,000 dilution

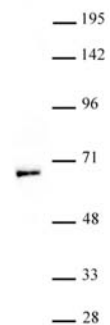
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at  $-20^{\circ}\text{C}$  for up to 2 years. Keep all reagents on ice when not 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.



**MOF / MYST1 antibody (pAb) tested by Immunofluorescence.**

Formaldehyde fixed HeLa cells stained with MOF / MYST1 antibody at a 1:2,000 dilution.



**MOF / MYST1 antibody (pAb) tested by Western blot.**

Whole-cell extract of HeLa cells (30  $\mu$ g) probed with MOF / MYST1 pAb at a dilution of 1:500.