

## Uhrf1 antibody (mAb)

**Catalog No:** 61341

**Clone:** 1RC-1C10

**Isotype:** IgG1, k

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

**Reactivity:** Human

**Quantity:** 100 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 0.93 µg/µl

**Molecular Weight:** 96 kDa

**Background:** Uhrf1 (ubiquitin-like, containing PHD and RING finger domains 1) is a transcription and cell cycle regulator. It functions as a putative E3 ubiquitin-protein ligase. Uhrf1 is important for G1/S transition of the cell cycle. It is also believed that Uhrf1 may help recruit the DNA methyltransferase, DNMT1, to hemimethylated DNA and assist in the stabilization of DNMT1 with chromatin.

**Immunogen:** This antibody was raised against a synthetic peptide corresponding to amino acid residue(s) 263-793 of human Uhrf1.

**Buffer:** Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 32 mM glycine, 0.07 mM EDTA, 30% glycerol, and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

Applications Validated by Active Motif:

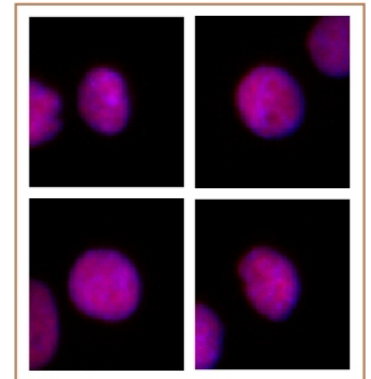
ICC/IF: 0.5 µg/ml dilution

WB\*: 1:1000 - 1:2,500 dilution

\*Note: Many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot. In addition, we recommend the addition of 0.05% Tween 20 to all blocking solutions to reduce background.

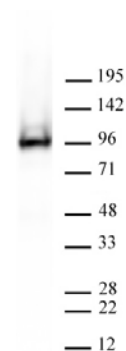
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This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



### Uhrf1 antibody (mAb) tested by immunofluorescence.

Formaldehyde fixed HeLa cells stained with Uhrf1 antibody at a 0.5 µg/ml dilution.



### Uhrf1 antibody (mAb) tested by Western blot.

20 µg of HeLa nuclear extract was run on SDS-PAGE and probed with antibody at 1:1000 dilution.