

## Ubn1 antibody (pAb)

**Catalog Nos:** 39869, 39870

**RRID:** AB\_2614993

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human

**Volumes:** 100 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 145 kDa

**Background:** UBN1 (Ubinuclein 1, also known as HIRA binding protein) is a ubiquitously expressed nuclear protein that acts as a novel regulator of senescence. UBN1 is involved in the formation of senescence-associated heterochromatin foci (SAHF), which represses expression of proliferation-promoting genes. It is expressed in a number of cancer cell lines and tumors and may be required for replication-independent chromatin assembly. UBN1 is part of a large multi-subunit protein complex that interacts with CEBPA and Epstein-Barr virus BLZF1. It co-localizes with HIRA, PML and SP100 in the PML bodies of senescent cells. When overexpressed, UBN1 accumulates in the nucleus of cells defective for cytokinesis.

**Immunogen:** This Ubinuclein antibody was raised against a recombinant protein corresponding to amino acids 1-345 of human Ubn1.

**Buffer:** Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 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:

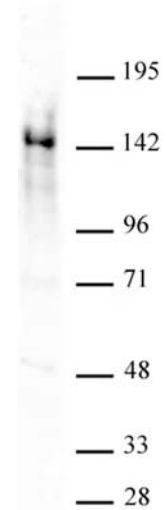
WB\*: 1:500 - 1:1,000 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

\*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.

**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.



**Ubn1 antibody (pAb) tested by Western blot.**  
Nuclear extract of HeLa cells (40 µg) probed with Ubn1 antibody at a dilution of 1:500.