

## Histone H1.0 antibody (pAb)

**Catalog Nos:** 61417, 61418

**RRID:** AB\_2793627

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human, Mouse

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

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 30 kDa

**Background:** Histone H1 is a linker histone, present at the interface between the nucleosome core and DNA entry/exit points. Histone H1 is responsible for establishing higher-order chromatin structure. **Histone H1.0** is a unique variant considered a replacement linker histone which is expressed and incorporated into chromatin in the absence of DNA replication. Histone H1.0 is more highly expressed in cells that are growth inhibited or terminally differentiated.

**Immunogen:** This antibody was raised against a peptide within the N-terminal region of human Histone H1.0.

**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:

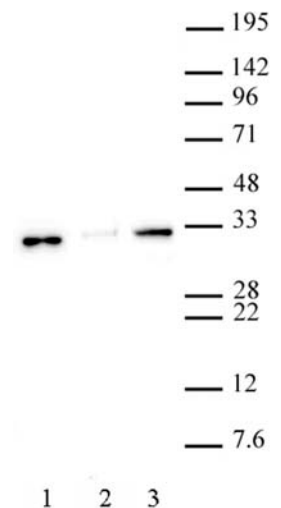
WB\*: 1:500 - 1:2,000 dilution

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.



### Histone H1.0 (pAb) tested by Western blot.

Nuclear Extract of 3T3-L1 cells (20  $\mu$ g) following stimulation with a defined hormonal mixture (DMSO, dexamethasone, insulin, and IBMX) were probed with Histone H1.0 (pAb) at a dilution of 1:500.

Lane 1: Recombinant human Histone H1.0 (100 ng).

Lane 2: Day 0 - preadipocytes.

Lane 3: Day 12 - fully differentiated adipocytes.