

## SUV420H1 antibody (pAb)

**Catalog No:** 61415, 61416

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human

**Purification:** Affinity Purified

**Host:** Rabbit

**Volume:** 100 µl, 10 µl

**Molecular Weight:** 100 kDa

**Background:** SUV420H1 (suppressor of variegation 4-20 homolog 1) is a histone methyltransferase that specifically trimethylates 'Lys-20' of histone H4. H4 'Lys-20' trimethylation represents a specific tag for epigenetic transcriptional repression. Mainly functions in pericentric heterochromatin regions, thereby playing a central role in the establishment of constitutive heterochromatin in these regions. SUV420H1 is targeted to histone H3 via its interaction with RB1 family proteins (RB1, RBL1 and RBL2) (By similarity).

**Immunogen:** This antibody was raised against a peptide within the N-terminal region of human SUV420H1.

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

**Application Notes:**

Validated Applications:

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.

Published Applications:

WB

**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Upon receipt, unconjugated antibodies may be stored at -20°C for up to 2 years. Fluorophore- & enzyme-conjugated antibodies should be stored at 4°C. Fluorophore-conjugated antibodies should be protected from light. Keep reagents on ice when not in storage; to avoid repeated freeze/thaw cycles, we recommend aliquoting items that will be stored frozen into single-use fractions prior to freezing. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

— 195

— 142

 — 96

— 71

— 48

**SUV420H1 antibody (pAb) tested by Western blot.**

Nuclear extract of Hep G2 cells (30 µg) probed with SUV420H1 antibody (pAb) at a dilution of 1:500.