

EuHMT1 antibody (pAb)

Catalog Nos: 39777, 39778

RRID: AB_2793339

Isotype: IgG

Application(s): WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Rabbit

Concentration: 1 µg/µl

Molecular Weight: 155 kDa

Background: EuHMT1 (Euchromatic Histone Methyltransferase 1, EHMT1, EuHMTase1, G9a like protein, GLP1) is a SET domain-containing histone methyltransferase specific for lysine 9 of histone H3. It is possible that this protein is involved in the silencing of Myc and E2F-responsive genes and thus could play a role in regulating cell cycle transitions from G0 to G1.

Immunogen: This EuHMT1 antibody was raised against a peptide derived from the N-terminal region of human EuHMT1.

Buffer: PBS pH 7.5 containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

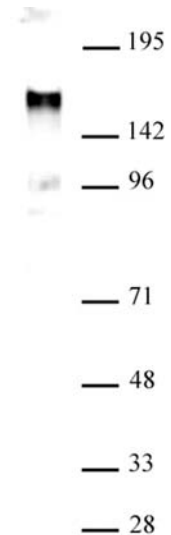
Applications Validated by Active Motif:

WB*: 1 - 2 µg/ml 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.

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.



EuHMT1 antibody (pAb) tested by Western blot.

Nuclear extract of HeLa cells (20 µg) probed with EuHMT1 antibody (1 µg/ml dilution).