

PHF20 antibody (pAb)

Catalog Nos: 61123, 61124

RRID: AB_2615013

Isotype: IgG

Application(s): WB

Reactivity: Human, Mouse

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 120 kDa

Background: PHF20 (PHD finger protein 20, Glioma-expressed antigen 2, TZP) is putative transcription factor and a component of the MLL complex. PHF20 contains a PHD finger, two Tudor domains, and interacts with histone H4 methylated at lysine 20.

Immunogen: This PHF20 antibody was raised against a peptide derived from human PHF20.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

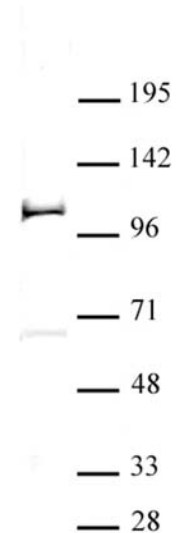
Application Notes:

Applications Validated by Active Motif:

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

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.



PHF20 antibody (pAb) tested by Western blot
20 μ g nuclear extract of F9 retinoic acid - treated mouse teratocarcinoma stem cells (Cat. No. 36113) probed with PHF20 antibody (1:500).