

PHF6 antibody (pAb)

Catalog Nos: 61807, 61808

RRID: AB_3216380

Application(s): WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Isotype: IgG

Molecular Weight: 45 kDa

Background: PHF6 is a member of the plant homeodomain finger (PHF) family, which has a potential role in regulation of chromatin accessibility to lineage-specific transcription factors as part of the nucleosome remodeling and deacetylase (NuRD) epigenetic complex. PHF6 has been reported as a tumor suppressor in association with hematopoietic malignancies including T-cell acute lymphoblastic leukemia (T-ALL) and acute myeloid leukemia (AML). PHF6 may have tissue-specific expression, and is overexpressed in both breast and colorectal cancers. Germline mutations in PHF6 are the cause of the Börjeson–Forssman–Lehmann X-linked intellectual disability syndrome (BFLS).

Immunogen: This antibody was generated using a peptide from the N-terminal region of human PHF6 (PHD finger protein 6 isoform 1).

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 - 1:1,000 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.