

PHF8 antibody (pAb)

Catalog Nos: 39711, 39712

RRID: AB_2793316

Isotype: Serum

Application(s): WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: None

Host: Rabbit

Molecular Weight: 135 kDa

Background: PHF8 (PHD finger protein 8) is a member of the JmjC-containing (Jumonji-C) class of histone demethylase proteins that are involved in the regulation of genome function through the removal of methyl groups from histones. PHF8 also has an N-terminal PHD finger and demethylates Histone H3 monomethyl Lys9 (H3K9me1) and Histone H3 dimethyl Lys9 (H3K9me2), Histone H3 dimethyl Lys27 (H3K27me2) (which are all modifications associated with transcriptional repression) and also Histone H3 dimethyl Lys36 (H3K36me2). Defects in PHF8 lead to Siderius type X-linked mental retardation (MRXSSD).

Immunogen: This PHF8 antibody was raised against a synthetic peptide derived from human PHF8.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

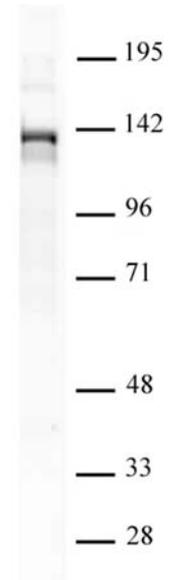
Application Notes:

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.



PHF8 pAb tested by Western blot.
Nuclear extract of HeLa cells (20 µg) was probed with PHF8 pAb (1:500 dilution).