

CBX8 antibody (pAb)

Catalog Nos: 61237, 61238

RRID: AB_2793563

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 45 kDa

Background: **CBX8** (chromobox homolog 8, PC3, RC1) is a Component of the Polycomb group (PcG) multiprotein PRC1 complex, a complex required to maintain the transcriptionally repressive state of many genes, including Hox genes, throughout development. PcG PRC1 complex acts via chromatin remodeling and modification of histones; it mediates monoubiquitination of histone H2A 'Lys-119', rendering chromatin heritably changed in its expressibility.

Immunogen: This CBX8 antibody was raised against a peptide corresponding to the C-terminal region of human CBX8 (Chromobox protein homolog 8).

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

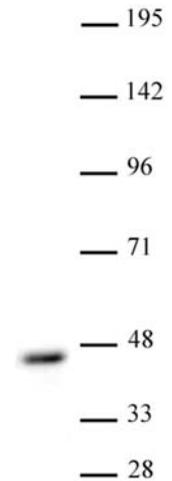
Application Notes:

Applications Validated by Active Motif:

WB: 1:500 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.



CBX8 antibody (pAb) tested by Western blot.

Whole-cell extract of HeLa cells (20 μ g) probed with CBX8 antibody (1:500).