

PBX1b antibody (mAb)

Catalog Nos: 61165, 61166

RRID: AB_2793534

Clone: CT 41.1

Isotype: IgG1, k

Application(s): ChIP, IHC, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 40 kDa

Background: PBX proteins are a family of TALE-class transcription factors that associate with a subset of Hox proteins, enhancing their DNA binding affinities and specificities. PBX proteins also heterodimerize with the Meis/Prep subfamily of to form trimeric complexes with Hox proteins at specific DNA sites. Amino acids 348-430 are absent in isoform **PBX1b**.

Immunogen: This PBX1b antibody was raised against a recombinant protein corresponding to amino acids 319 - 347 of human PBX1b.

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

Application Notes:

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.