

## PBX2 antibody (mAb)

---

**Catalog Nos:** 61155, 61156

**RRID:** AB\_2793530

**Clone:** 2.1

**Isotype:** IgG1

**Application(s):** IHC, WB

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 51 kDa

---

**Background:** PBX proteins are a family of TALE-class transcription factors that 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. PBX2 is the predominant high-molecular-weight isoform PBX protein detectable by immunoblotting and is widely expressed during embryonic development, particularly in neural and epithelial tissues.

**Immunogen:** This PBX2 antibody was raised against a recombinant protein corresponding to amino acids 308 - 430 of mouse PBX2.

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

**Application Notes:**

This antibody is also available as an AbFlex® engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91227.

**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.