

PITX2 antibody (pAb)

Catalog Nos: 61303, 61304

RRID: AB_2614945

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 33 kDa

Background: PITX2 (paired-like homeodomain transcription factor 2, ARP1, RGS, RIEG, RIEG1) controls cell proliferation in a tissue-specific manner and is involved in morphogenesis. During embryonic development, exerts a role in the expansion of muscle progenitors. May play a role in the proper localization of asymmetric organs such as the heart and stomach.

Immunogen: This PITX2 antibody was raised against a peptide within the N-terminal region of human PITX2.

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

Application Notes:

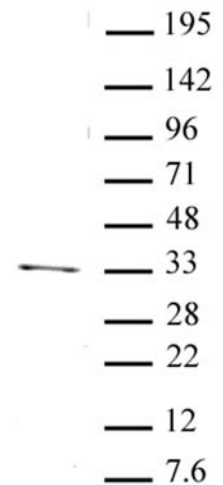
Validated Applications:

WB: 1:500 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

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.



PITX2 antibody (pAb) tested by Western blot.
Whole cell extract of Saos-2 cells (20 μ g) probed with PITX2 antibody (1:500).