

Prox1 antibody (mAb)

Catalog Nos: 61091, 61092

RRID: AB_2615008

Clone: 4G10

Isotype: IgG2b

Application(s): IF, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 90 kDa

Background: Prox1 (Prospero-like homeobox protein 1) is a sequence-specific DNA binding transcription factor. Prox1 controls lymphatic endothelial and retinal progenitor cell differentiation, as well as development of the liver and lens of the eye. Required for lymphatic endothelial cells specification and the formation of lymphatic vessels.

Immunogen: This Prox1 antibody was raised against a recombinant protein corresponding to amino acids 547-737 of human Prox1.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 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.