

PAX9 antibody (mAb)

Catalog Nos: 61077, 61078

RRID: AB_2793498

Clone: 7C2

Isotype: IgG1

Application(s): IHC, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 38 kDa

Background: **PAX9** (PAX-9, Paired box protein 9) is a sequence specific DNA binding transcription factor required for normal development of the thymus, parathyroid glands, teeth, portions of the skull and larynx as well as distal limbs. **PAX9** is a member of the paired box (PAX) family of transcription factors, which typically contain a paired box domain, a paired type homeodomain and an octapeptide sequence.

Immunogen: This PAX9 antibody was raised against a recombinant protein corresponding to amino acids 263-342 of mouse PAX9 protein.

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.