

PAX7 antibody (mAb)

Catalog Nos: 39803, 39804

RRID: AB_2793348

Clone: 2F12H4

Isotype: IgG2b

Application(s): ICC, IF, WB

Reactivity: Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 55 kDa

Background: PAX7 (Paired-box protein 7, PAX7) is a member of the paired box (PAX) family of transcription factors. These proteins are DNA binding transcriptional activators and typically contain a paired box domain, an octapeptide, and a paired-type homeodomain. These genes play critical roles during mammalian development and cancer growth. The specific function of the PAX7 gene is unknown but speculated to play a role in myogenesis. PAX7 is involved in the specification of satellite cells in skeletal muscle. PAX7 can bind to DNA as a heterodimer with Pax-3.

Immunogen: This PAX7 antibody was raised against a peptide derived from mouse PAX7.

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.