

Myf-5 antibody (mAb)

Catalog Nos: 39801, 39802

RRID: AB_2614936

Clone: 1E2G4

Isotype: IgG2a

Application(s): ICC, IF, WB

Reactivity: Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 28 kDa

Background: Myf-5 (Myogenic Factor 5, MFY5) is a basic helix-loop-helix transcription factor involved in muscle differentiation. Myf-5 induces differentiation of fibroblasts into myoblasts. Efficient DNA binding requires dimerization with another bHLH protein.

Immunogen: This Myf-5 antibody was raised against a peptide derived from mouse Myf-5.

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.