

## MyoD antibody (mAb)

---

**Catalog Nos:** 39991, 39992

**RRID:** AB\_2793421

**Clone:** 5F11

**Isotype:** IgG2a

**Application(s):** ICC, IF, WB

**Reactivity:** Human, Mouse

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Rat

**Concentration:** 1 µg/µl

**Molecular Weight:** 40 kDa

---

**Background:** MyoD (Myogenic differentiation factor 1, Myf-3), is a transcription factor belonging to the myogenic factors family that contains a basic helix-loop-helix (bHLH) DNA-binding domain. MyoD is the central regulator in the signaling pathways that regulate muscle development.

**Immunogen:** This MyoD antibody was raised against a recombinant protein corresponding to full-length mouse MyoD.

**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.