# Myosin (human) monoclonal antibody (F36.5B9)

This antibody is covered by our Worry-Free Guarantee.

Citations: 10

View Online »

## **Ordering Information**

Order Online »

ALX-BC-1040-S-L001	1ml
ALX-BC-1040-S-L005	5ml

Manuals, SDS & CofA

View Online »

### **Handling & Storage**

**Use/Stability** Stable for a few weeks when stored at +4°C.

**Handling** Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

#### **Product Details**

Application IHC (FS), WB

Clone F36.5B9

Formulation Liquid. In sterile filtered supernatant with 0.09% sodium

azide.

**Host** Mouse

**Immunogen** Human ventricular myosin.

lgG2a

**Recommendation Dilutions/Conditions**Immunohistochemistry (frozen sections, can be used

undiluted)Western Blot (dilution 1:10)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined

individually for each application.

Species Reactivity Human

**Specificity** Recognizes the S1/S2 junction of human ventricular

myosin heavy chain β.

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

-

Last modified: May 29, 2024



info-

eu@enzolifesciences.com