Myosin monoclonal antibody (F109.3E1)

This antibody is covered by our Worry-Free Guarantee.

Citations: 20

View Online »

Ordering Information

Order Online »

ALX-BC-1150-S-L001	1ml
ALX-BC-1150-S-L005	5ml

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC (FS), WB

Clone F109.3E1

Formulation Sterile filtered supernatant containing 0.1% sodium azide.

Host Mouse

Immunogen Both human ventricular myosin light chains.

lsotype lgG1

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, Rat

Specificity Recognizes myosin light chain 2 (LC2) (aa 45-59).

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

.

Last modified: May 29, 2024

