

Muscle actin monoclonal antibody (HHF35) (Ready-to- Use)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-30931	~6 µg - 6ml
-----------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC (PS), WB

Clone HHF35

Formulation Liquid. *Ready-to-use* solution in PBS containing 0.1% sodium azide and 1% BSA.

Host Mouse

Immunogen Human myocardium (proteins extracted in 2% SDS).

Isotype IgG1

Purity Detail Affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections; this antibody can be used for staining of paraffin-embedded tissue sections fixed in neutral buffered formalin and is compatible with most commonly used histological fixatives. Choose a good quality enzyme-linked anti-mouse IgG1 antibody for detection)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Monkey, Rabbit

Specificity Recognizes actin isotypes α and γ of skeletal, cardiac and smooth muscle. This antibody labels rhabdomyosarcomas, hemangiopericytomas, leiomyosarcomas and leiomyomas but does not react with non-myogenic sarcomas. It is nonreactive with other neoplasms such as carcinoma or lymphoma and with other mesenchymal and all epithelial cells, with the exception of myoepithelium.

UniProt ID P68133 (ACTA1), P63267 (ACTG2)

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)