

# $\beta$ -Actin polyclonal antibody

Actin is one of the major proteins of the cytoskeleton. It regulates contractile potential in muscle tissues and helps to control the shape as well as the motility of non-muscle cells. Actin is expressed in all eukaryotic cells. In higher eukaryotes, there are six isoforms of actin.

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

Citations: 7

[View Online »](#)

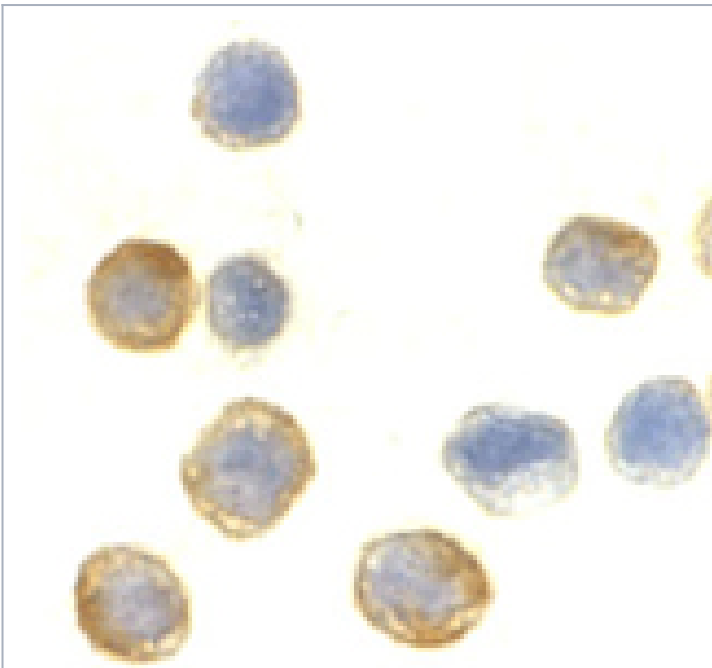
## Ordering Information

[Order Online »](#)

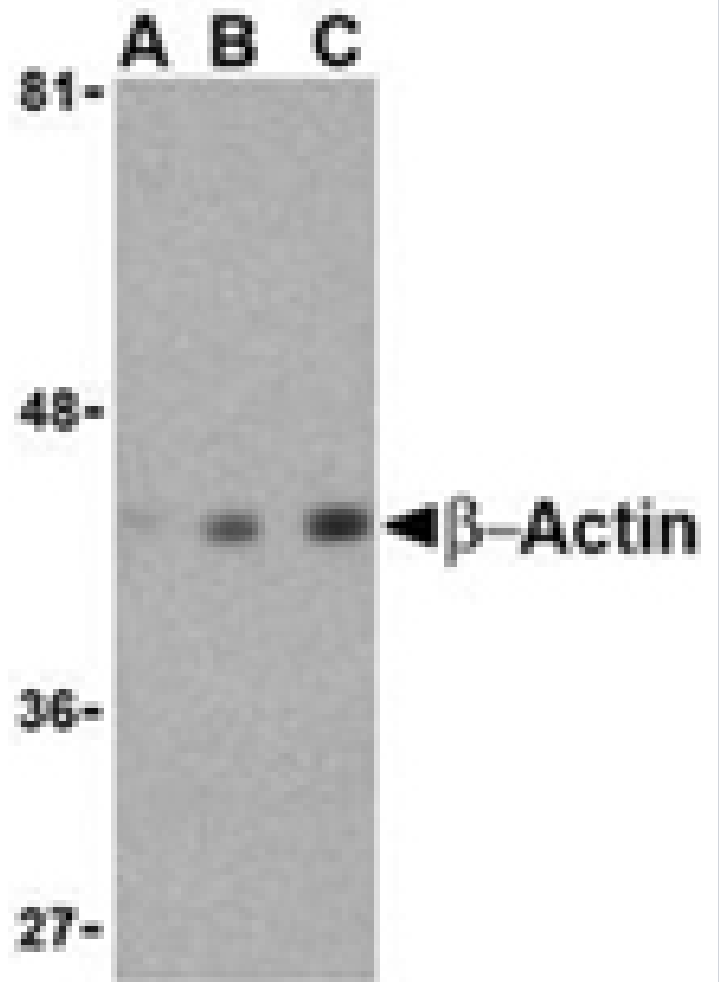
ADI-905-733-100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Immunocytochemistry analysis of beta-actin in HeLa cells with beta-actin antibody at 10 $\mu$ g/ml.



Western blot analysis of beta-actin in HeLa cell lysate with beta-actin antibody at (A) 0.5, (B) 1 and (C) 2 $\mu$ g/ml.

## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	ACTA, ACTSA
Application	ICC, WB
Application Notes	Detects a band of ~43kDa by Western blot.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the sequence near the C-terminus of human $\beta$ -Actin.
Purity Detail	Peptide affinity purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (10 $\mu$ g/ml)Western Blot (1-2 $\mu$ g/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	P60709
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



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-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

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