

# Titin monoclonal antibody (F146.9B9)

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

Citations: 3

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-BC-3010-S-L001	1ml
ALX-BC-3010-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

Alternative Name Connectin

Application IHC (FS), WB

Clone F146.9B9

Formulation 1ml/5ml of sterile filtered supernatant containing 0.1% sodium azide.

Host Mouse

Immunogen Purified titin from human skeletal muscle.

Isotype IgG1

**Recommendation** Immunohistochemistry (frozen sections)Western Blot (dilution 1:10)  
**Dilutions/Conditions** Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse, Rabbit

UniProt ID Q8WZ42

**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)