

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 Dilutions/Conditions | Immunohistochemistry (frozen sections)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, Mouse, Rabbit |
| UniProt ID | Q8WZ42 |
| Worry-free Guarantee | This antibody is covered by our Worry-Free Guarantee |

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)