

# Ryanodine receptor monoclonal antibody (34C)

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

Citations: 8

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-804-016-R100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** RyR

**Application** ICC, IHC (FS), IP, WB

**Application Notes** Detects a band of ~565kDa by Western blot, representing RyR from rat skeletal muscle extracts. In non-mammalian vertebrates, a doublet is seen at 565kDa, representing the  $\alpha$  and  $\beta$  isoforms.

**Clone** 34C

**Formulation** Liquid. Contains 0.05% sodium azide.

**Host** Mouse

**Immunogen** Partially purified chicken pectoral muscle RyR (ryanodine receptor).

**Isotype** IgG1

**Recommendation Dilutions/Conditions** Immunohistochemistry (frozen sections, 1:1,000)Western Blot (1:5,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Amphibian, Bovine, Dog, Fish, Human, Mink, Monkey, Mouse, Rabbit, Rat, Sheep

**Specificity** Recognizes RyR-1 and RyR-2 isoforms. In chicken, it detects the  $\alpha$ ,  $\beta$  and cardiac isoforms; in frog the  $\alpha$  and  $\beta$  isoforms; and in fish the  $\alpha$  isoform. Also recognizes RyR-3 in mouse cells.

UniProt ID

Q90984

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-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)