

Ryanodine receptor monoclonal antibody (C3-33)

The ryanodine receptor (RyR) is the channel responsible for the release of calcium from the sarcoplasmic reticulum (SR) in muscle cells and also plays a role in calcium regulation in non-muscle cells. The RyR exists as a homotetramer and is predicted to have a short cytoplasmic C-terminus and 4-10 transmembrane domains; the remainder of the protein, termed the "foot" region is located in the cytoplasm between the T-tubule and the SR. The mammalian RyR is the product of three different genes: RyR-1, which is expressed predominantly in skeletal muscle and areas of the brain, RyR-2, which is expressed predominantly in the heart muscle but also found in the stomach, endothelial cells and diffuse areas of the brain, and RyR-3 which is found in smooth muscle and the brain (striatum, thalamus and hippocampus).

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

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|------------------|-------|
| ALX-804-015-C100 | 100µg |
|------------------|-------|

Manuals, SDS & CofA

[View Online »](#)

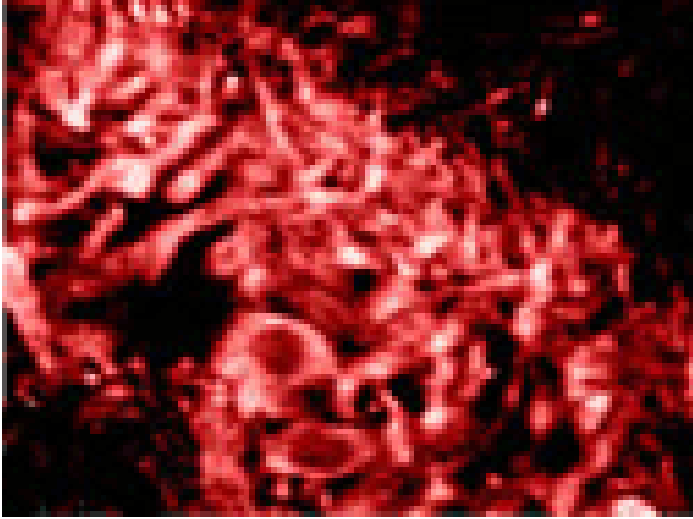


Figure: Immunolocalization of ryanodine receptor 1 (RyR1) in rat brain hippocampus using Monoclonal Antibody to Ryanodine Receptor (C3-33) (Prod. No. ALX-804-015).

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 | Flow Cytometry, ICC, IHC (FS), IHC (PS), IP, WB |
| Application Notes | Detects a band of 565kDa by Western blot. In non-mammalian vertebrates, a doublet is seen at 565kDa representing the α and β isoforms of the receptor. |
| Clone | C3-33 |
| Crossreactivity | Cross-reacts weakly with RyR-1. |
| Formulation | Liquid. In PBS containing 0.05% sodium azide. |
| Host | Mouse |
| Immunogen | Dog cardiac RyR (ryanodine receptor). |
| Isotype | IgG1 |
| Recommendation | Immunocytochemistry (1:100-1:200)Immunohistochemistry (frozen sections, |
| Dilutions/Conditions | 1 μ g/ml)Immunohistochemistry (paraffin sections, 1 μ g/ml)Western Blot (1 μ g/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application. |
| Species Reactivity | Amphibian, Chicken, Dog, Fish, Guinea pig, Human, Mouse, Rabbit, Rat |
| UniProt ID | Q92736 (human RYR2) |



ENZO LIFE SCIENCES
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com
be@enzolifesciences.com

France
Phone: +33 1 72 44 01 555
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

This antibody is covered by our Worry-Free Guarantee