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 »

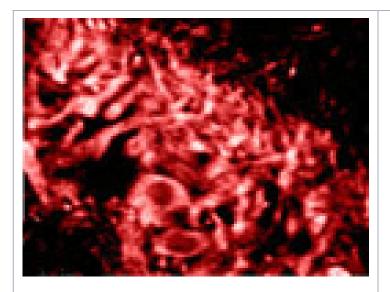


Figure: Immunolocalization of ryanodine receptor 1 (RyR1) in rat brain hippocampus using Monoclonal Antibody to Ryanodine Receptor (C3-33) (Prod. No. ALX-804-015).

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

Cross-reacts weakly with RyR-1.

Formulation Liquid. In PBS containing 0.05% sodium azide.

Host Mouse

Immunogen Dog cardiac RyR (ryanodine receptor).

lsotype lgG1

Recommendation Dilutions/Conditions Immunocytochemistry (1:100-1:200)Immunohistochemistry

(frozen sections, 1µg/ml)Immunohistochemistry (paraffin sections, 1µg/ml)Western Blot (1µg/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)

Last modified: May 29, 2024



info-

eu@enzolifesciences.com