

# Inositol 1,4,5- trisphosphate receptor (type-I) polyclonal antibody

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

## Ordering Information

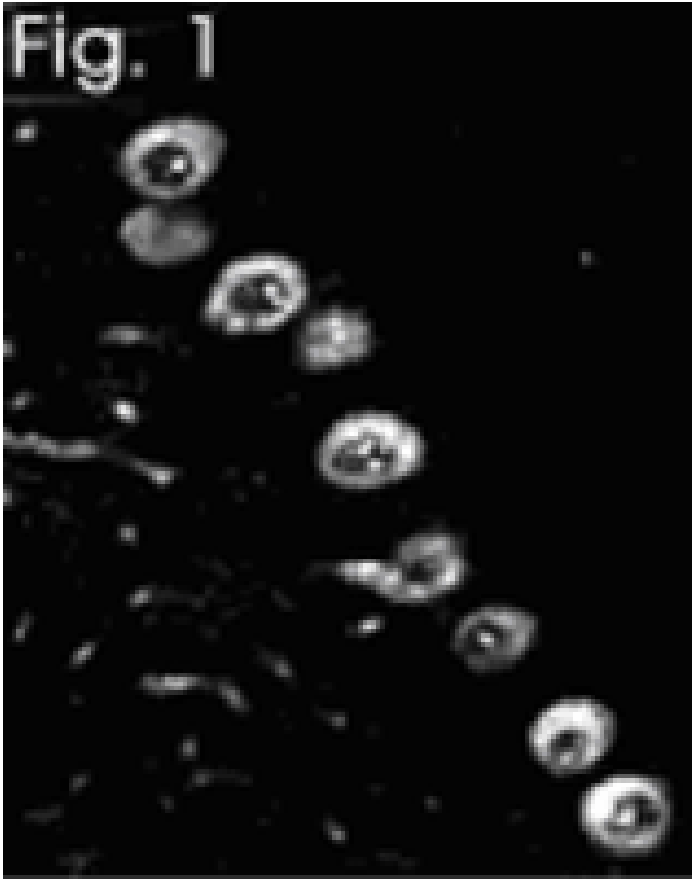
[Order Online »](#)

ALX-210-169-R100	100µl
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)

Fig. 1



**Figure:** Immunolocalization of IP<sub>3</sub>R-I in rat cerebellum using PAb to Inositol 1,4,5-Trisphosphate Receptor (Type 1)

## 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** IP3 receptor type-I, IP3R-I, ITPR1, INSP3R1, IP3R1

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

**Application Notes** Detects a band of ~240kDa by Western blot.

**Crossreactivity** Weakly cross-reacts with human IP3R-I.

**Formulation** Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to aa 1829-1848 N1829KKKDDEVDRDAPSRKKAKE1848) of the cytoplasmic C-terminal domain of human IP3R-I (inositol 1,4,5-triphosphate receptor (type-I)). This sequence is completely conserved in mouse and rat IP3R-I.

**Purity Detail** Epitope-affinity purified IgG.

**Recommendation Dilutions/Conditions** Immunohistochemistry (frozen sections, 1:250)Western Blot (1:1,000)Optimal conditions must be determined individually for each application.

**Species Reactivity** Dog, Mouse, Rat

**UniProt ID** Q14643

**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)