

# Calreticulin polyclonal antibody

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

Citations: 4

[View Online »](#)

## Ordering Information

[Order Online »](#)

|                  |       |
|------------------|-------|
| ALX-210-126-R100 | 100µl |
|------------------|-------|

Manuals, SDS & CofA

[View Online »](#)

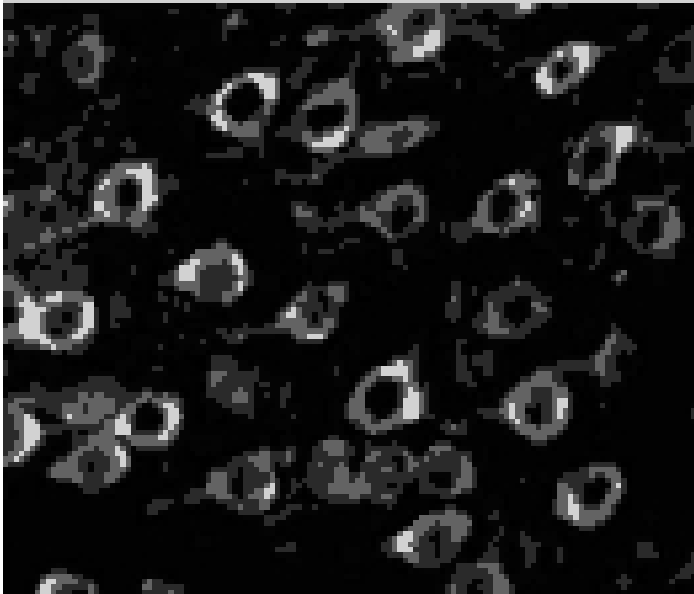


Figure: Immunolocalization of calreticulin in rat brain cortex using PAb to calreticulin (Prod. No. [ALX-210-126](#)).

Calreticulin polyclonal antibody IHC, Calreticulin polyclonal antibody Immunohistochemistry, yes

ALX-210-127, 210-

127.gif, [http://static.enzolifesciences.com/fileadmin/files/image/210-](http://static.enzolifesciences.com/fileadmin/files/image/210-127.gif)

127.gif, **Figure:** Immunocytochemistry of  $\beta$ -COP in NIH-3T3 cells using PAb to  $\beta$ -COP (Prod. No. [ALX-210-127](#)).

).

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** Flow Cytometry, ICC, IHC, IP, WB

**Formulation** Liquid. Antiserum containing 0.05% sodium azide.

**Host** Rabbit

**Immunogen** Recombinant human calreticulin.

**Recommendation Dilutions/Conditions** Immunohistochemistry (1:100)Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Dog, Human, Monkey, Mouse, Rabbit, Rat

**UniProt ID** P27797

**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)