

# Calreticulin (NT) polyclonal antibody

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

## Ordering Information

[Order Online »](#)

ALX-210-170-R100	100µl
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)

~66 kDa



**Figure:** Western blot of calreticulin in AtT-20 cell lysates using PAb to Calreticulin (NT) (Prod. No. ALX-210-170).

## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	CALR, CRT
Application	ICC, WB
Application Notes	Detects a band of ~60-66kDa by Western blot.
Formulation	Liquid. Contains 1mg/ml BSA and 0.05% sodium azide.
Host	Chicken
Immunogen	Synthetic peptide corresponding to aa 24-43 (K24EQFLDGDWNTNRWVESKHK43) of mouse calreticulin.
Purity Detail	Epitope-affinity purified IgY.
Recommendation Dilutions/Conditions	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, Hamster, Human, Mouse, Rat
UniProt ID	P14211
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

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-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@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)