

Calreticulin polyclonal antibody

Calreticulin (CRT) is a multifunctional, multi-compartmental protein most abundant in the ER lumen. CRT contains the ER-retrieval sequence, KDEL, and has been best characterized as a soluble molecular chaperone of new or misfolded proteins and a Ca^{2+} -binding protein. Both CRT and its membrane bound homolog, Calnexin (CNX) interact with proteins and glycoproteins that have monoglucosylated N-glycans. The CRT/CNX cycle promotes correct folding, inhibits aggregation of folding intermediates, blocks premature oligomerization, regulates ER degradation, and provides quality control by preventing incompletely folded glycoproteins from exiting to the Golgi complex.

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

Citations: 36

[View Online »](#)

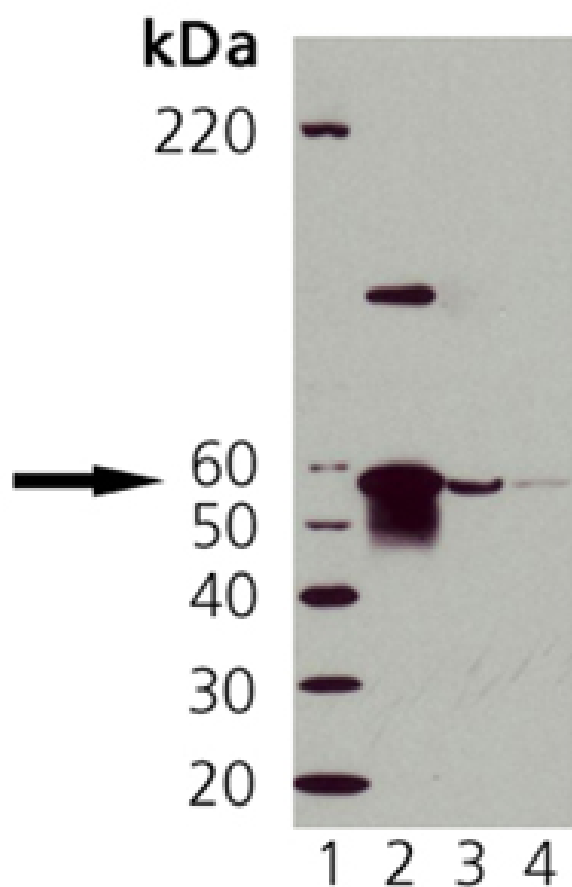
Ordering Information

[Order Online »](#)

ADI-SPA-600-J	1mg
ADI-SPA-600-D	50µg
ADI-SPA-600-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western Blot Analysis of Calreticulin: Lane 1: MWM,
Lane 2: HeLa cell lysate, Lane 3: PC-12 cell lysate.
Lane 4: 3T3 cell lysate

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	CALR, CRT
Application	ICC, IHC, IP, WB
Application Notes	Predicted MW of ~63kDa.
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
GenBank ID	AD000092
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the sequence near the C-terminus of human Calreticulin.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000, human; 1:500, mouse; ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Bovine, Dog, Human, Mouse, Rat
UniProt ID	P27797
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)