Calreticulin polyclonal antibody

The multifunctional, multi-compartmental protein Calreticulin (Crt) functions as a soluble molecular chaperone of new or misfolded proteins as well as a Ca2+-binding protein. Most abundant in the ER lumen, Crt expression also occurs in other membrane-bound organelles, the cell surface, and extracellularly. Also known as CRP-55, calregulin and HACBP (high affinity calcium-binding protein), Crt contains the ER-retrieval sequence, KDEL, and is the soluble paralog of the ER membrane protein Calnexin (Cnx). Crts three domains include a 180 residue N-terminal domain, a proline-rich P-domain (residues 189-288) that binds Ca2+ with high affinity and shares homology with Cnx and calmegin, and a 110 residue C-terminal domain that binds Ca2+ with low affinity but high capacity. The P-domain may interact with the co-chaperone ERp57 (Grp58), a thiol reductase. The NMR structure of the P-domain consists of an extended hairpin that appears to form a curved protrusion from the Crt core domain. Both Crt and its membrane bound homolog CNX interact with proteins and glycoproteins possessing monoglucosylated N-glycans. The Crt/Cnx cycle promotes correct folding, inhibits aggregation of folding intermediates, blocks premature oligomerization, regulates ER degradation, and prevents incompletely folded glycoproteins from exiting to the Golgi complex. Crt also appears to function as an auto-antigen in systemic lupus erythematosus, rheumatoid arthritis, celiac disease, complete congenital heart block, and halothane hepatitis. A diversity of additional functions attributed to Crt includes adhesion, blood function, and cardiac and neuronal development gene expression.

This antibody is covered by our Worry-Free Guarantee.

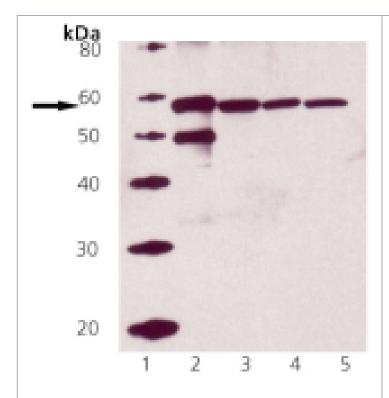
Ordering Information

Order Online »

ADI-SPA-602-D	50µg
ADI-SPA-602-F	200μg

Manuals, SDS & CofA

View Online »



Western blot analysis of Calreticulin pAb: Lane 1: MW marker, Lane 2: 50 ng Calreticulin, Lane 3: HeLa, Lane 4: 3T3, Lane 5: PC12.

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 WB

Application Notes Detects a band of ~55kDa by Western blot.

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium

azide.

Rabbit Host

Immunogen Recombinant human calreticulin.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:1,000, 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 Human, Mouse, Rat

UniProt ID P27797

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 eu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 fr@enzolifesciences.com

Phone: +49 7621 5500 526 de@enzolifesciences.com

Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 uk@enzolifesciences.com