

# Calnexin (luminal domain) (human), (recombinant)

Calnexin (CNX), an abundant ~90 kDa molecular chaperone, is a resident type I transmembrane protein of the endoplasmic reticulum (ER) (ref.1,2). A majority of the calnexin protein resides in the luminal portion of the ER, and is the membrane-bound paralog of the soluble ER protein Calreticulin. In mammalian cells, calnexin and calreticulin (CRT) are lectins that play key roles in glycoprotein folding within the ER, specifically binding oligosaccharide intermediates that contain a single terminal GlcManGlcNAc residue. Calnexin associates with newly synthesized, incompletely folded monomeric glycoproteins as well as numerous oligomeric protein complexes, including

Citations: 1

[View Online »](#)

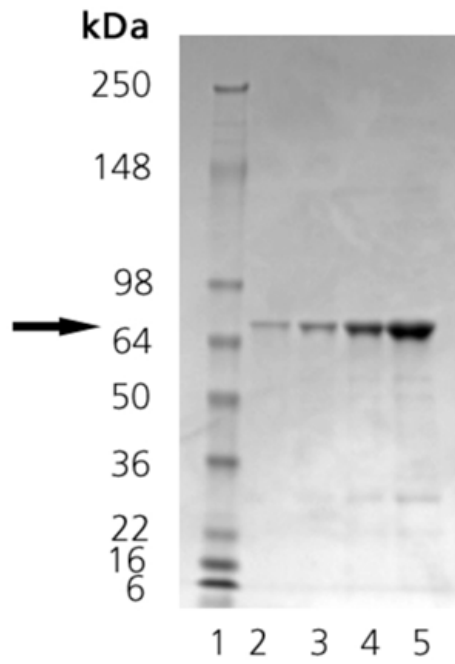
## Ordering Information

[Order Online »](#)

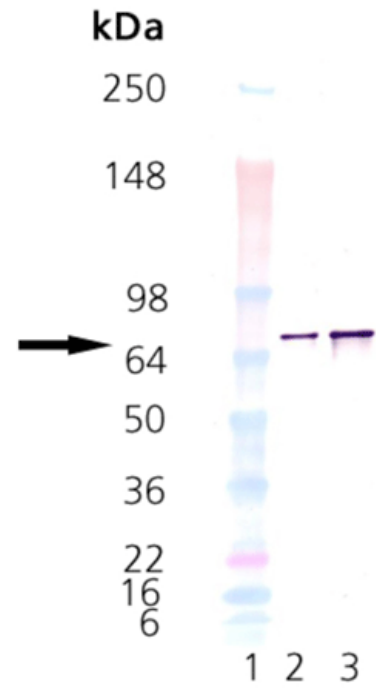
ADI-SPP-865-D	50µg
---------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis of Calnexin luminal domain (Prod. No. ADI-SPP-865) stained with Coomassie. Lane 1: MWM, Lane 2: 0.5 $\mu$ g; Lane 3: 1 $\mu$ g; Lane 4: 2 $\mu$ g; Lane 5: 5 $\mu$ g



Western Blot analysis of Calnexin luminal domain (Prod. No. ADI-SPP-865) probed with 1:1,000 dilution of Prod. No. ADI-SPA-865. Lane 1: MW marker, Lane 2: 50ng; Lane 3: 100ng

## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	CNX
Application Notes	Western blot control.
Formulation	Liquid. In TRIS containing sodium chloride, 5% glycerol, and 5mM calcium chloride.
MW	~74kDa
Purity	≥85% (SDS-PAGE; Western blot) (full length protein)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> .
UniProt ID	P27824

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)