

# Calnexin monoclonal antibody (AF18)

Calnexin, also referred to as IP90 or p88, is a ~88kDa chaperone protein found in the endoplasmic reticulum. It binds with partially folded immunoglobulin proteins and retains them in the endoplasmic reticulum until they are completely folded.

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

Citations: 16

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## 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** p88, IP90

**Application** ICC, IP, WB

**Application Notes** Detects a band of ~90kDa by Western blot.

**Clone** AF18

**Formulation** Liquid. Ascites diluted in PBS containing 0.05% sodium azide.

**Host** Mouse

**Immunogen** Focus human hepatoma cell line.

**Isotype** IgG1

**Recommendation Dilutions/Conditions** Immunocytochemistry (1:500) Immunoprecipitation, Western Blot (1:4,000) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Species Reactivity** Human, Mouse

**UniProt ID** P27824

**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)

Last modified: May 29, 2024