

# Cpn10 polyclonal antibody

Chaperonin 10 (Cpn10 or Hsp10), is the ~10 kDa mammalian equivalent to GroES of E. coli. Chaperonins play fundamental roles in the folding, assembly, and translocation of other proteins. Chaperonin 10 has been located in human platelets but more importantly it may also be present in human maternal serum. It has been reported that human Cpn10 is identical with Early pregnancy factor (EPF), which is involved in control over cell growth and development.

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

Citations: 10

[View Online »](#)

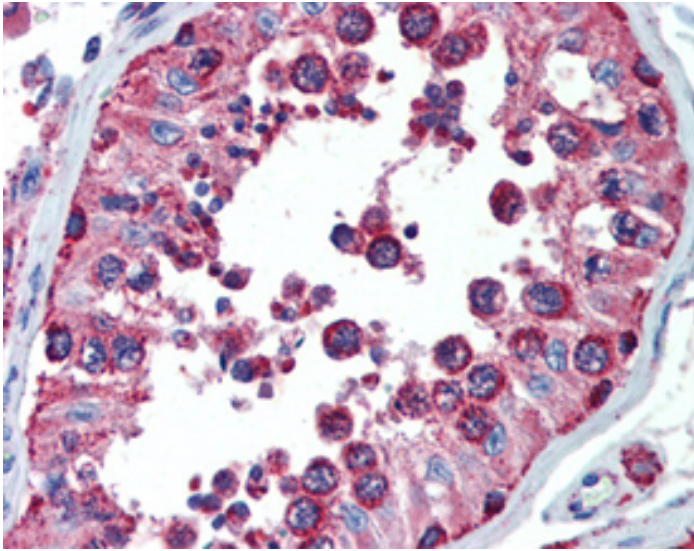
## Ordering Information

[Order Online »](#)

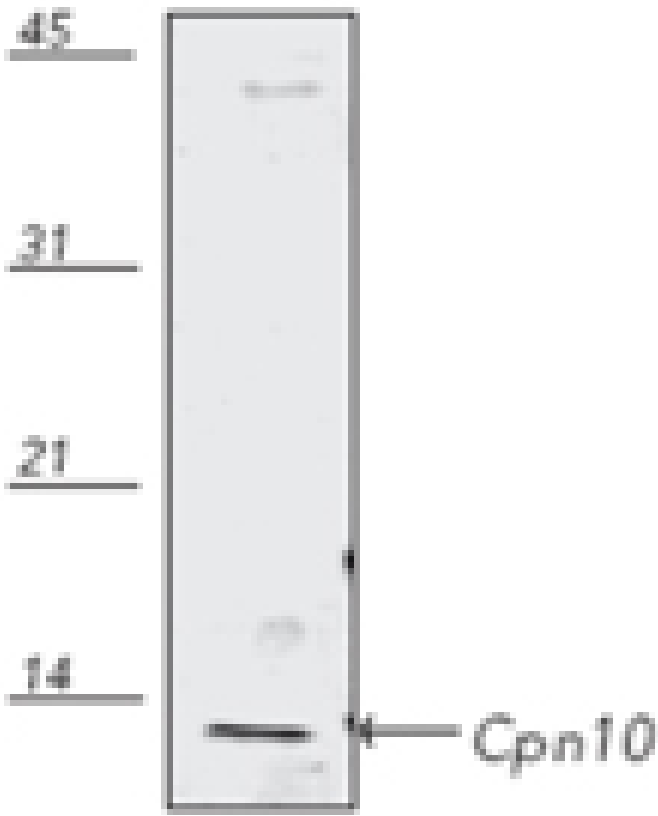
ADI-SPA-110-D	50µg
ADI-SPA-110-F	200µg

## Manuals, SDS & CofA

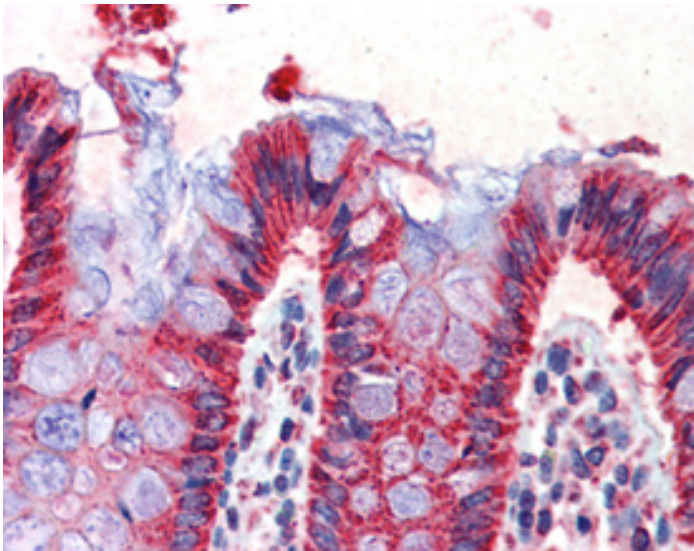
[View Online »](#)



IHC Analysis: Human testis tissue stained with Cpn10, pAb at 10µg/ml.



Western blot analysis of human HeLa (heat shocked) cell lysate, probed with Cpn10 (Hsp10) pAb.



IHC Analysis: Human colon tissue stained with Cpn10, pAb at 10µg/ml.

# 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	HSP10, HSPE1, GroES, Heat shock protein 10, Chaperonin 10
Application	IHC (PS), WB
Application Notes	Detects a band of ~10kDa by Western blot.
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
Gene/Protein Identifier	NM_002157 (RefSeq)
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the sequence near the C-terminus of human Cpn10.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:2,000, 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	Bovine, Dog, Guinea pig, Human, Mouse, Porcine, Rabbit, Rat, Sheep, Xenopus
Technical Info / Product Notes	Recommended by the Human Protein Atlas Organization for IHC (Ensembl No. ENSG00000115541).
UniProt ID	P61604

Worry-free Guarantee

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

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)