

# Cpn10 (human), (recombinant)

Chaperonin 10 (Cpn 10) is the ~10kD mammalian equivalent to GroES of *E. coli*. Chaperonins play fundamental roles in the folding, assembly, and translocation of other proteins. Chaperonin function in vivo requires two distinct family members, Cpn60 and Cpn10, both of which are also present in chloroplasts and mitochondria. The best studied Cpn60 and Cpn10, are the GroEL and GroES proteins of *E. coli*, respectively. Cpn10 and Cpn60 both exhibit a sevenfold axis of symmetry and function as a team in the protein folding and assembly processes. 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 to early pregnancy factor (EPF), which is involved in control over cell growth and development. This identification suggests that Cpn10 may act like a hormone in stressful situations such as pregnancy.

Citations: 2

[View Online »](#)

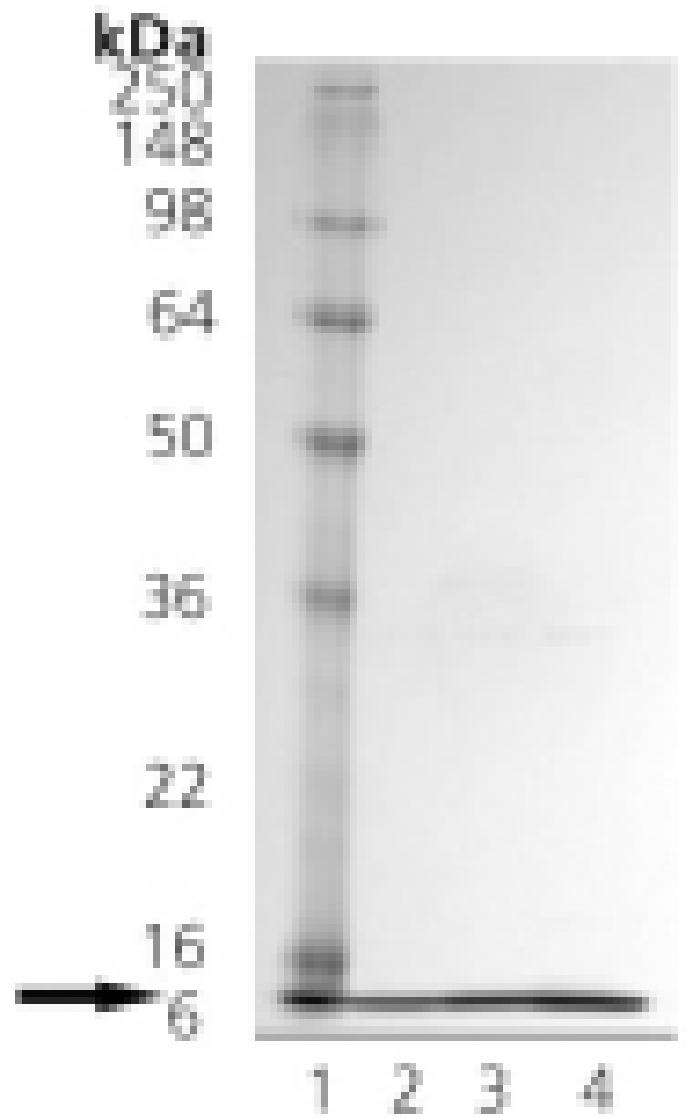
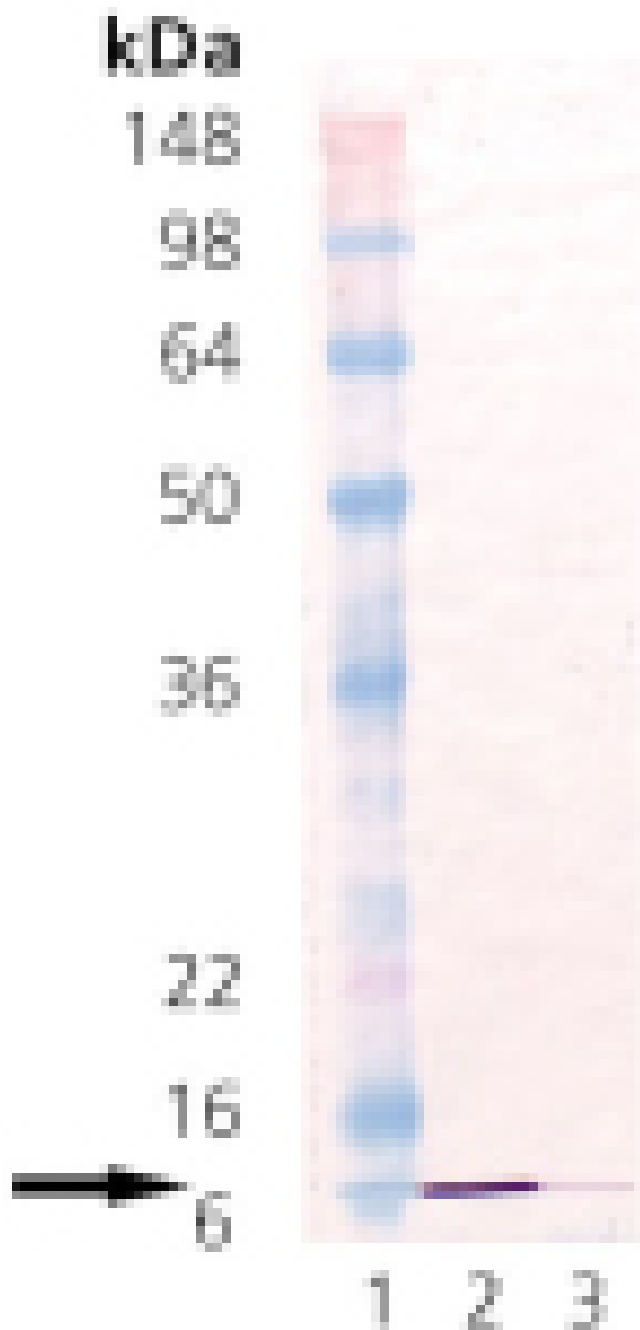
## Ordering Information

[Order Online »](#)

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

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage      -80°C

Shipping      Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** HSP10, HSPE1, GROES, Heat shock protein 10, Chaperonin 10

**Application Notes** Western blot control.

**Formulation** Liquid. In 20mM TRIS, pH 8.0, containing 150mM sodium chloride.

**MW** ~10kDa

**Purity** ≥90% (SDS-PAGE; Western blot)

**Purity Detail** Purified by multi-step chromatography.

**Source** Produced in *E. coli*.

**UniProt ID** P61604

**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)