

Cpn10 (human), (recombinant)

Recombinant human Cpn10 that assists protein folding and assembly as a mitochondrial and cytosolic chaperonin.

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 »](#)

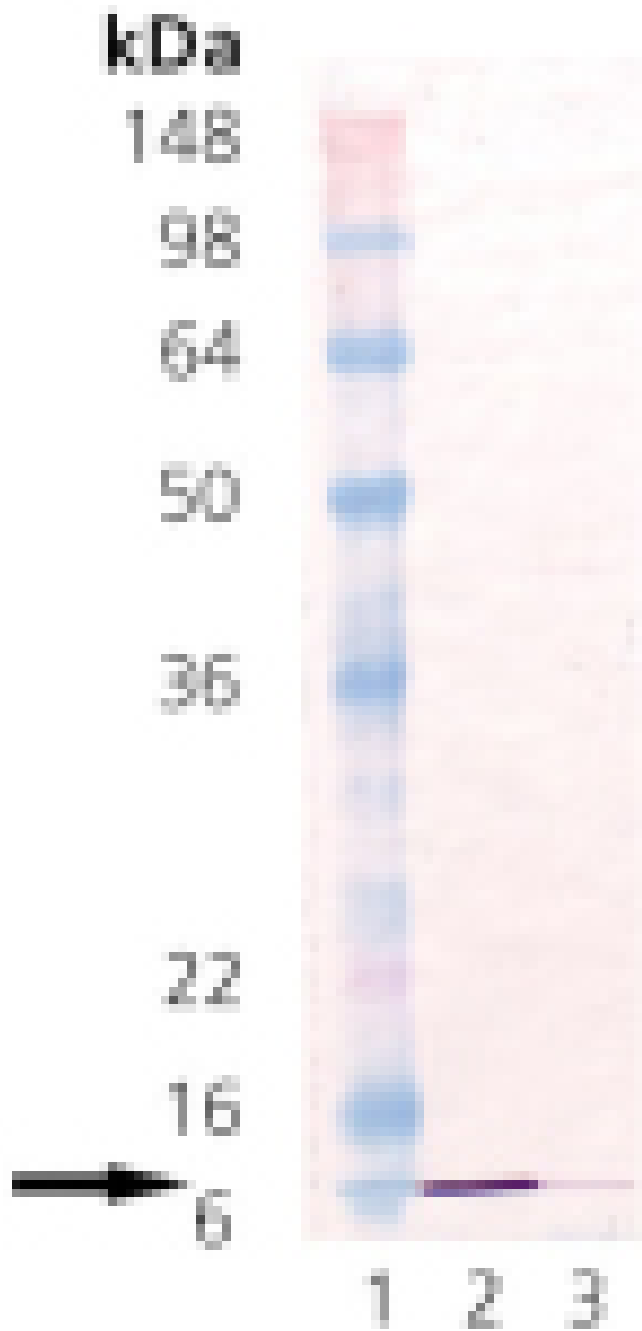
Ordering Information

[Order Online »](#)

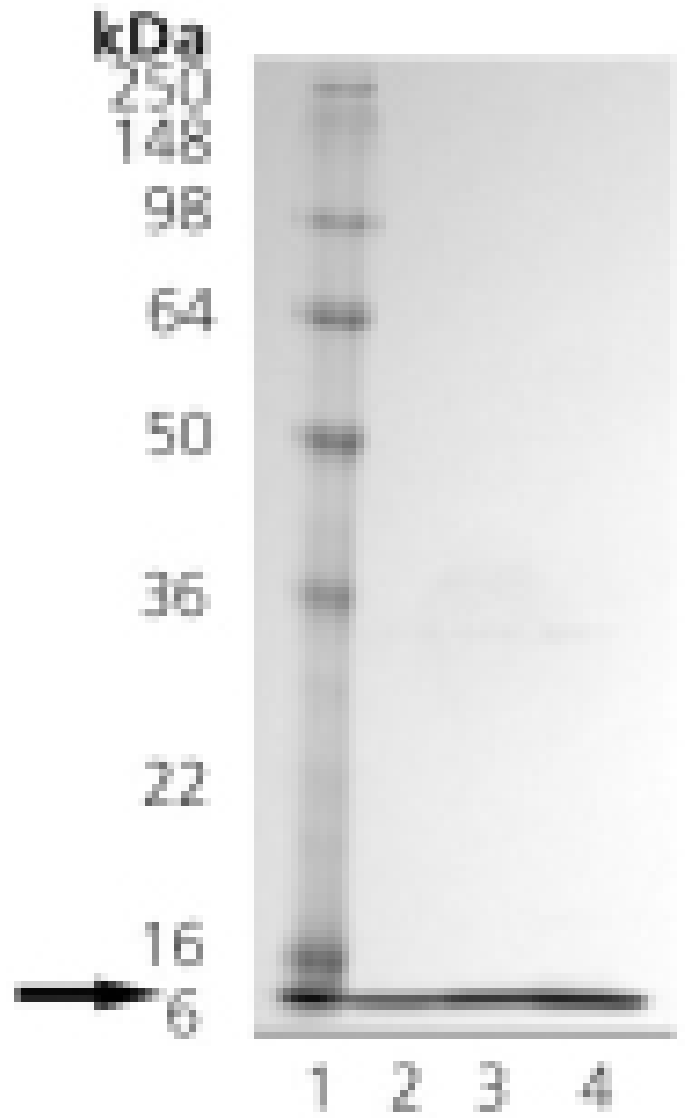
ADI-SPP-110-F	200µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis: Lane 1: MW marker, Lane 2: Cpn10 Recombinant Human Protein, Lane 3: HeLa lysate (Prod. No. ADI-LYC-HL100) probed with Cpn10 polyclonal antibody (Prod. No. ADI-SPA-110).



SDS-PAGE analysis: Lane 1: MWM, Lane 2: 1 μ g, Lane 3: 2 μ g, Lane 4: 5 μ g Cpn10 Recombinant Human Protein.

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 <i>E. coli</i> .
UniProt ID	P61604

Last modified: October 9, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

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

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

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