## 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 »

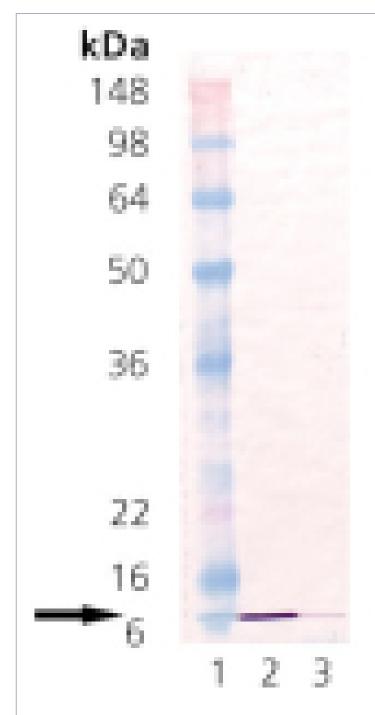
## **Ordering Information**

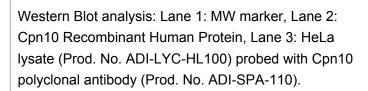
**Order Online** »

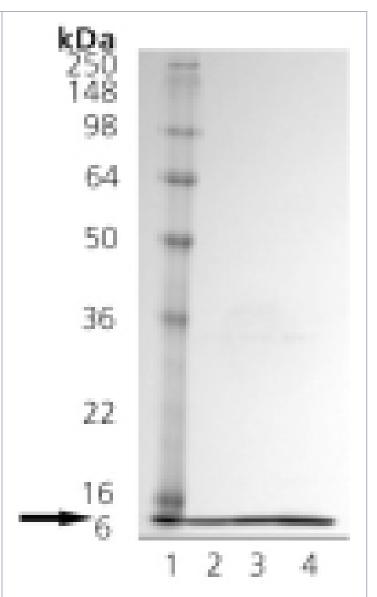
ADI-SPP-110-D	50µg
ADI-SPP-110-F	200µg

Manuals, SDS & CofA

View Online »







SDS-PAGE analysis: Lane 1: MWM, Lane 2: 1µg, Lane 3: 2µg, Lane 4: 5µg 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 *E. coli*.

UniProt ID P61604

Last modified: October 9, 2025

