HOP (human), (recombinant)

Hsp70-Hsp90 Organizing Protein (HOP, p60) is an ~60kDa protein that is a critical intermediate component for the efficient maturation of steroid receptor complexes, serving to recruit Hsp90 to Hsp70-containing complexes. Unactivated steroid hormone receptors are found in heterooligomeric complexes that are thought to stabilize a partially folded receptor polypeptide prior to hormone - dependent activation. Hsp70 and Hsp90 are both known to associate with steroid hormone receptors. Numerous studies have shown that Hsp70 functions in an ATPdependent manner through transient interactions to mediate folding or unfolding of polypeptide chains. Hsp90 is thought to perhaps also function in some capacity related to folding or protein-protein interactions. HOP contains three tetratricopeptide repeat (TPR) domains, TPR1, TPR2a and TPR2b. Hsp70 binding has been localized to TPR1, and sp90 binding has been localized to TPR2a. Importantly, the highly conserved EEVD sequence that terminates many Hsp70 family members, and the similar MEEVD sequence that terminates Hsp90, are important recognition sites for the TPR domains. The co-crystal structures for TPR1 plus a GPTIEEVD octapeptide (Hsp70 sequence) and TPR2a plus the MEEVD pentapeptide (Hsp90) have been solved. HOP is closely related to a human 63 kDa protein that is sensitive to simian virus SV40 transformation and is related to the yeast heat shock-responsive STI1 gene product.

Citations: 1

View Online »

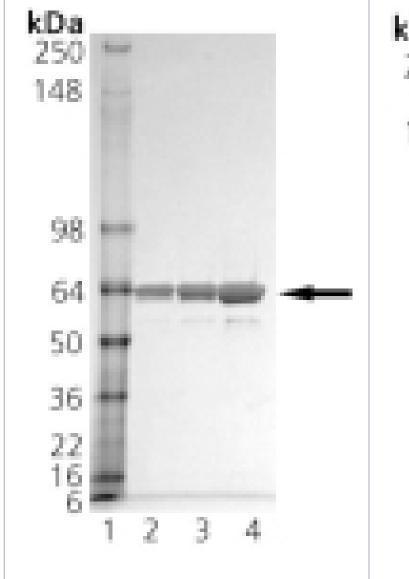
Ordering Information

Order Online »

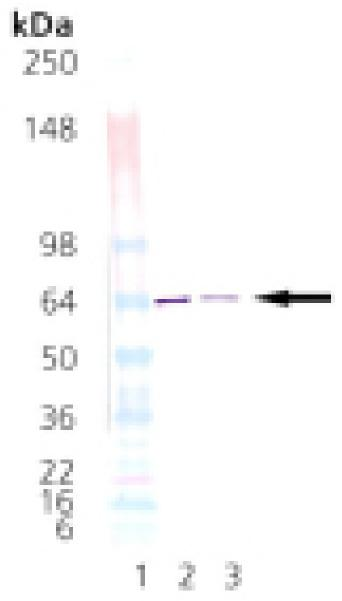
ADI-SRP-1510-D	50µg
ADI-SRP-1510-F	200µg

Manuals, SDS & CofA

View Online »



SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 1µg, Lane 3: 2µg, Lane 4: 5µg HOP Recombinant Protein.



Western Blot analysis: Lane 1: MWM, Lane 2: HOP Recombinant Protein, Lane 3: HeLa cell lysate.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name HSC70/HSP90-organizing protein, STIP1, STI1, Stress-

induced phosphoprotein 1, STUB1, p60, CHIP

Application Notes Western blot control.

Formulation Liquid. In Dulbecco's PBS.

MW ∼63kDa

Purity ≥90% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in *E. coli*.

UniProt ID P31948

Last modified: May 29, 2024



eu@enzolifesciences.com