HSP60 (mouse), (recombinant)

Recombinant mouse Hsp60, a mitochondrial chaperone that assists in protein folding and oligomer assembly, induced by stress and linked to immune and inflammatory

Mouse Hsp@Sbelongs to a highly conserved family that includes molecular chaperones from a variety of species such as plant Hsp60 (known as Rubisco binding protein), GroEL, the E. coli Hsp60, and the 65 kDa major antigen of mycobacteria. Hsp60 is localized in the mitochondrial matrix of eukaryotes, and in the chloroplast of plants. Hsp60s from divergent species share a number of common characteristics: high abundance; induction with environmental stress such as heat shock; homo - oligomeric structures of either 7 or 14 subunits which reversibly dissociate in the presence of Mg and ATP; ATPase activity; and a role in folding and assembly of oligomeric protein structures. These similarities correspond with studies in which the single -ring human mitochondrial homolog Hsp60 and its co - chaperonin Hsp10 were expressed in an E. coli strain engineered to keep the groE operon under strict regulatory control. The findings demonstrate that expression of Hsp60 -Hsp10 enabled successful performance of all essential in vivo functions of GroEL and its co-chaperonin, GroES. Several studies reveal a possible link between members of the Hsp60 family and a number of autoimmune diseases, atherosclerosis, and chlamydial disease. Overexpression of self Hsp60 is seen in the synovial tissue of rheumatoid arthritic (RA) patients, and can accompany both cellular and humoral reactivity against Hsp60 in RA. Chlamydial heat shock protein Hsp60, a homolog of E. coli GroEL, appears capable of eliciting macrophage activation, and several studies reveal a correlation between Hsp60 responses and the immunopathologic manifestations of human chlamydial disease.

Citations: 3

View Online »

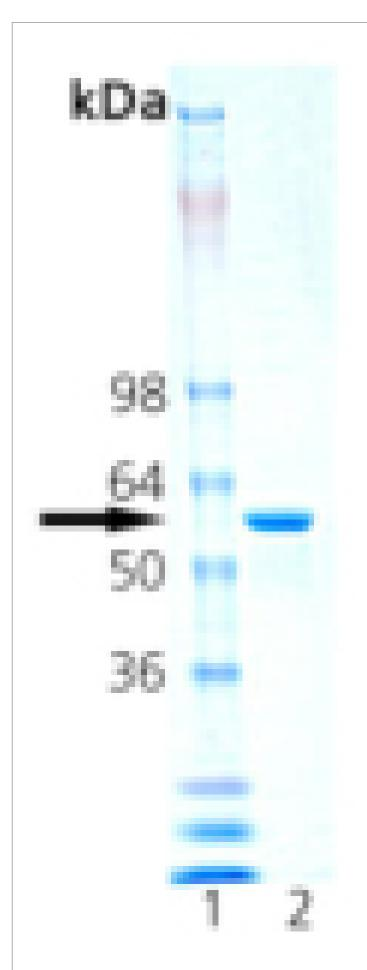
Ordering Information

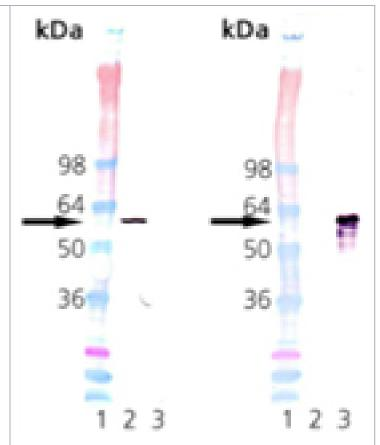
Order Online »

ADI-SPP-741-D	50µg
ADI-SPP-741-F	200μg

Manuals, SDS & CofA

View Online »





Western Blot analysis: Lane 1: MW marker; Lane 2: 100ng Mouse HSP60 Protein (Prod. No. ADI-SPP-741); Lane 3: 100ng E. coli GroEL Protein (Prod. No. ADI-SPP-610); Left Blot probed with HSP60 mAb (Prod. No. ADI-SPA-806) at 1.0μg/ml; Right Blot probed with GroEL mAb (Prod. No. ADI-SPS-870) at 1.0μg/ml.

SDS-PAGE analysis: Lane 1: MWM, Lane 2: 2.0µg of purified Mouse HSP60 Protein (Prod. No. ADI-SPP-741)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Chaperonin 60, CPN60, HspD1, Heat shock protein 60

Application NotesATPase activity assay (positive). Western blot control.

Formulation Liquid. In 10mM TRIS-HCl, pH 7.5, containing 150mM

sodium chloride, 1.0mM EDTA, and 1.0mM DTT.

MW ~60kDa

Purity ≥90% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in *E. coli*.

UniProt ID P63038

Last modified: October 9, 2025

