

HSP60 Protein

Full-length recombinant protein expressed in *E. coli* cells

Catalog # H33-54G

Lot # C211-2

Product Description

Recombinant full-length human HSP60 was expressed in *E. coli* cells using an N-terminal GST tag. The gene accession number is [NM_002156](#).

Gene Aliases

HSPD1, CPN60, GROEL, HSP65, SPG13, HuCHA60

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

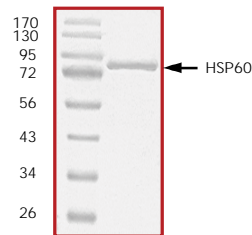
HSP60 is a chaperonin protein required for ATP-dependent folding of precursor polypeptides and complex assembly. HSP60 prevents aggregation and mediates protein refolding after heat shock. Mutations and variations in the HSP60 gene product may be implicated in genetic diseases (1). HSP60 has been strongly implicated as an example of molecular mimicry in the pathogenicity of autoimmune diseases in T cell-mediated protection. Human HSP60 was found to activate the complement system in normal human serum in a dose-dependent manner. Since complement activation plays an important role in the development of atherosclerosis, the levels of complement-activating anti-HSP60 antibodies are elevated in atherosclerosis-related diseases. (2)

References

1. Pochon, N A. et al: Genetic complexity of the human hsp 60 gene. *Int. Immunol.* 1996; 8(2):221-30
2. Zoltán, P. et al: Antibodies against human heat-shock protein (hsp) 60 and mycobacterial hsp65 differ in their antigen specificity and complement-activating ability. *Int. Immunol.* 1999;11 (9), 1363-1370.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

Purity



The purity was determined to be **>95%** by densitometry.
Approx. MW **85kDa**.

HSP60 Protein

Full-length recombinant protein expressed in *E. coli* cells

Catalog Number H33-54G

Specific Lot Number C211-2

Purity	>95%
Concentration	0.2 $\mu\text{g}/\mu\text{l}$
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.