

HSP10 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog # H29-54G

Lot # G022-1

Product Description

Recombinant full-length human HSP10 was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM_002157](#).

Gene Aliases

CPN10; GROES; HSPE1

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

HSP10 is a 10-kDa stress-inducible mitochondrial protein which is a member of the heat shock proteins. HSP10 is synthesized at elevated rates in cultured rat hepatoma cells challenged with heat shock or amino acid analogues. HSP10 has been implicated in multiple disease states in humans. In myocardial samples from patients with chronic atrial fibrillation, cardiomyocytes respond to chronic atrial fibrillation with increased expression of HSP10 and HSP60 (1). HSP10 is also overexpressed early in prostate carcinogenesis. HSP10 is also overexpressed in large bowel carcinomas and it may have diagnostic and prognostic significance in the management of this tumour (2).

References

1. Schafner, A E. et al: Overexpression of heat shock protein 60/10 in myocardium of patients with chronic atrial fibrillation. *Ann Thorac Surg.* 2002 Sep;74(3):767-70.
2. Cappello, F. et al: The expression of HSP60 and HSP10 in large bowel carcinomas with lymph node metastase. *BMC Cancer.* 2005 Oct 28;5:139.

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 **>90%** by densitometry.
Approx. MW **36kDa**.

HSP10 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog Number H29-54G

Specific Lot Number G022-1

Purity	>90%
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.