

## H11 (HSP22) Protein

Full-length recombinant human protein expressed in Sf9 cells

**Catalog # H09-34G**

Lot # A1147-3

### Product Description

Full-length recombinant human H11 (HSP22) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [NM\\_014365](#).

### Gene Aliases

HSP22; CRYAC; CMT2L; DHMN2; E2IG1; HSPB8; HMN2; HMN2A; PP1629

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 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.

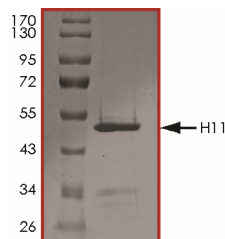
### Scientific Background

H11 or heat shock 22kDa protein 8 belongs to the superfamily of small heat-shock proteins containing a conservative alpha-crystallin domain at the C-terminal part of the molecule. The expression of H11 is induced by estrogen in estrogen receptor-positive breast cancer cells, and H11 also functions as a chaperone in association with Bag3, a stimulator of macroautophagy. H11 is involved in regulation of cell proliferation, apoptosis, and carcinogenesis (1). H11 also causes motor neuron-specific neurite degeneration (2).

### References

1. Benndorf, R.et.al: Shocking degeneration. Nature Genet. 36: 547-548, 2004.
2. Irobi, J.et.al: Mutant HSPB8 causes motor neuron-specific neurite degeneration. Hum. Molec. Genet. 19: 3254-3265, 2010.

### Purity



The purity of H11 (HSP22) protein was determined to be **>95%** by densitometry.  
Approx. MW **50 kDa**.

## H11 (HSP22) Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog #	H09-34G
Lot #	A1147-3
Purity	>95%
Concentration	0.1 µg/µl
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.

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

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**