HSP70B' (human), (recombinant)

Recombinant human HSP70B', an inducible 70 kDa stress protein that assists in protein folding, trafficking, and prevents aggregation under cellular stress conditions.

In response to adverse changes in their environment, cells from many organisms increase the expression of a class of proteins referred to as heat shock or stress proteins. One class of stress proteins is the HSP70 family which is comprised of a group of highly conserved, ubiquitous proteins of ~70 kDa that are located within different intracellular compartments and involved in a variety of cellular processes such as protein trafficking, folding, and prevention of aggregation. One strictly inducible member of the HSP70 family is Hsp70B'. In humans, the hsp70B' gene is 77% similar to hsp70 and it encodes a ~70 kDa protein which is more basic than Hsp70 (1). Hsp70 and Hsp70B' differ in their extent and pattern of inducibility in response to conditions such as heat shock and CdCl2 treatment (1). In humans, Hsp70 is constitutively expressed under normal conditions, but Hsp70B' is only induced in response to stress (1). There is no basal expression of Hsp70B' (1).

Citations: 2

View Online »

Ordering Information

Order Online »

ADI-SPP-762-E

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Hsp70B prime, Heat shock protein 70 B'

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

Formulation Liquid. In Dulbecco's PBS, pH 7.2, containing 8.1mM

sodium phosphate . 7H₂O, 1.5mM potassium phosphate, 2.7mM potassium chloride, and 137mM sodium chloride.

MW ~70kDa

Purity ≥85% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in *E. coli*.

Technical Info / Product NotesUS Patent No. 7,326,574.

UniProt ID P17066

Last modified: October 9, 2025

