

**HSP90b (HSP84)**

**CATALOG NO.:** HSP-21-619

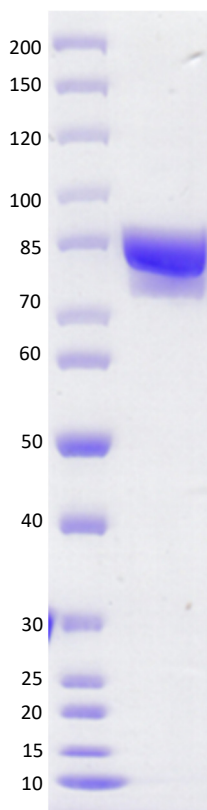
**LOT NO.:**

**DESCRIPTION:** Human recombinant HSP90b (residues 2-724; Genbank Accession # NM\_007355.4; MW = 85.65 kDa) expressed with an N-terminal StrepII tag an insect cell/baculovirus expression system.

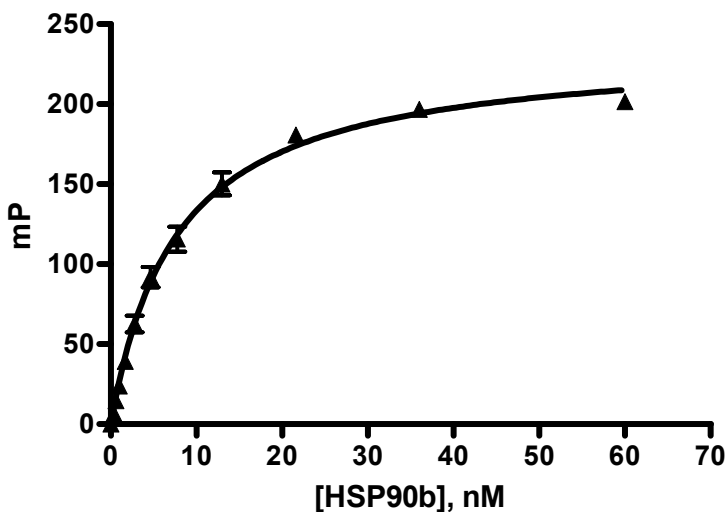
**PURITY:** >90% by SDS-PAGE.

**SUPPLIED AS:** \_\_\_ µg/µL in 50mM Tris pH8.0, 200mM NaCl, 10% glycerol, 1mM TCEP as determined by OD<sub>280</sub>.

**STORAGE:** -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice/ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.



**Coomassie blue-stained SDS-PAGE (8-16% acrylamide) of 4 µg of RBC HSP90b.** MW markers (left) are, from top, 200, 150, 120, 100, 85, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



**HSP90b Binding Assay.** FITC-Geldanamycin probe binding was detected using fluorescence polarization. The 20µL reaction contained 5nM probe and variable concentration of HSP90a. Polarization values (mP) were measured at an excitation wavelength at 485 nm and an emission wavelength at 520 nm using an PHERAstar reader (BMG Labtech) following a 3hr incubation.

This product is not intended for therapeutic or diagnostic use in animals or in humans.

**Reaction Biology**

1 Great Valley Parkway, Malvern PA, USA 19355  
 requests@reactionbiology.com www.reactionbiology.com