

StressXpress®  
HSP90 alpha  
ELISA Kit

93 /100 11 Citations

Powered by Bioz

[See more details on](#)



Discovery through Partnership | Excellence through Quality

Colorimetric detection of HSP90 Alpha  
Catalog No. SKT-107

### Product Name

HSP90 alpha ELISA Kit

### Description

Colorimetric detection of HSP90 Alpha

### Species Reactivity

Avian (Zebra finch), Goat, Human

### Platform

Microplate

### Sample Types

Cell lysates, Serum, Tissue, Whole Blood

### Detection Method

Colorimetric Assay

### Assay Type

Sandwich ELISA (Enzyme-linked Immunosorbent Assay)

### Utility

ELISA kit used to quantitate HSP90 alpha concentration in samples.

### Sensitivity

0.117 ng/ml

### Assay Range

0.44 - 28 ng/ml

### Incubation Time

30 minutes

## Number Of Samples

40 samples in duplicate

## Other Resources

Kit Booklet Lot No. SH531948 , Kit Booklet Lot No. SH187452 , Kit Booklet Lot No. SH388280 , Kit Booklet Lot No. KH387128 , MSDS

## Field Of Use

Not for use in humans. Not for use in diagnostics or therapeutics. For in vitro research use only.

## Properties

### Storage Temperature

4°C and -20°C

### Shipping Temperature

Blue Ice

### Product Type

ELISA Kits

### Assay Overview

1. Prepare Standard and samples in Standard and Sample Diluent. 2. Add 100 µL of Standard or sample to appropriate wells. 3. Cover plate with Plate Sealer and incubate at 37°C for 1 hour. 4. Wash plate four times with 1X Wash Buffer. 5. Add 100 µL of Biotinylated Antibody Working Solution to each well. 6. Cover plate with Plate Sealer and incubate at room temperature, 20-25 °C for 1 hour. 7. Wash plate four times with 1X Wash Buffer. 8. Add 100 µL of Streptavidin-HRP Working Solution to each well. 9. Cover plate with Plate Sealer and incubate at room temperature for 30 minutes. 10. Wash plate four times with 1X Wash Buffer. 11. Add 100 µL of TMB Substrate to each well. 12. Develop the plate in the dark at room temperature for 30 minutes. 13. Stop reaction by adding 100 µL of Stop Solution to each well. 14. Measure absorbance on a plate reader at 450 nm.

### Kit Overview

#### Component No.

#### Item

#### Quantity / Size

#### SKC-107A

Anti-Hsp90a Immunoassay Plate

1 Plate

**SKC-107B**

5X Hsp90a Extraction Reagent

1 vial/10 ml

**SKC-107C**

Recombinant Hsp90a Standard

2 vials

**SKC-107D**

Standard and Sample Diluent

1 vial/ 50 ml

**SKC-107E**

10X Wash Buffer Concentrate

1 vial/100 ml

**SKC-107F**

Anti-Hsp90a Biotinylated Antibody Concentrate

1 vial/150 µl

**SKC-107G**

Anti-Hsp90a Biotinylated Antibody Diluent

1 vial/ 13 ml

**SKC-107H**

Streptavidin: HRP Concentrate

1 vial/150 µl

**SKC-107I**

Streptavidin: HRP Diluent

1 vial/ 13 ml

**SKC-107J**

---

TMB Substrate

---

1 vial/ 13 ml

---

### **SKC-107K**

---

Stop Solution

---

1 vial/ 13 ml

---

### **Cite This Product**

---

HSP90 alpha ELISA Kit (StressMarq Biosciences Inc., Victoria BC CANADA, Catalog # SKT-107)

---

## **Biological Description**

---

### **Alternative Names**

HSP86, HSP89A, HSP90A, HSP90AA1, HSP90alpha, HSPC1, HSPCA, HsoCAL3

---

### **Research Areas**

---

Cancer, Cell Signaling, Chaperone Proteins, Heat Shock, Protein Trafficking, Tumor Biomarkers

---

### **Scientific Background**

---

Heat Shock Protein 90 (HSP90) is a highly conserved molecular chaperone essential for maintaining protein homeostasis across all eukaryotic cells. Although classified as a stress protein, HSP90 is abundantly expressed even under non-stress conditions, comprising up to 2% of cytosolic protein. Its core functions include the folding, maturation, stabilization, and trafficking of a wide range of client proteins.

HSP90 plays a pivotal role in regulating key signaling molecules such as kinases (v-Src, Wee1, c-Raf), transcription factors (p53), steroid receptors, and enzymes like telomerase and viral polymerases. These interactions are mediated through ATP-dependent conformational changes and co-chaperone complexes involving Cdc37, p23, and immunophilin-like proteins, which protect client proteins from proteasomal degradation.

In neuroscience, HSP90 is increasingly recognized for its role in modulating protein misfolding and aggregation—central features of neurodegenerative diseases such as Alzheimer's, Parkinson's, and Huntington's. Alterations in HSP90 expression or function can disrupt neuronal signaling and impair the activity of neuroprotective proteins, contributing to disease progression.

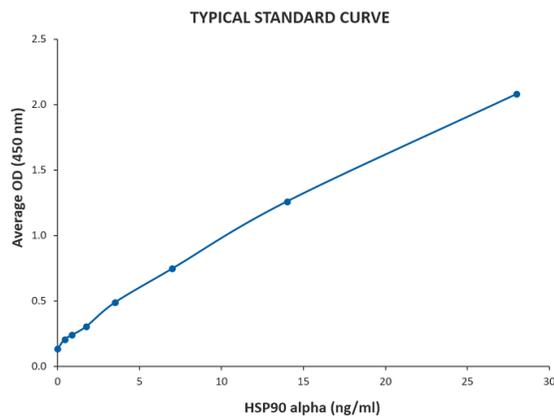
Pharmacological inhibitors of HSP90, including geldanamycin and radicicol, are being explored as therapeutic agents to restore proteostasis and reduce toxic protein accumulation in neurodegenerative models.

---

### **References**

1. Arlander S.J.H., et al. (2003) J Biol Chem. 278: 52572-52577.
2. Pearl H., et al. (2001) Adv Protein Chem. 59: 157-186.
3. Neckers L, et al. (2002) Trends Mol Med. 8:S55-S61.
4. Pratt W., Toft D. (2003) Exp Biol Med. 228:111-133.
5. Pratt W., Toft D. (1997) Endocr Rev. 18: 306-360.
6. Pratt W.B. (1998) Proc Soc Exptl Biol Med. 217: 420-434.
7. Whitesell L., et al. (1994) Proc Natl Acad Sci USA. 91: 8324- 8328.

## Product Images



Typical Standard Curve for the HSP90 Alpha ELISA kit (Enzyme-Linked Immunosorbent Assay) StressXpress® – SKT-107. Assay Type: Sandwich ELISA. Detection Method: Colorimetric Assay. Assay Range: 0.44 – 28 ng/mL.

## Reviews

Based on validation through cited publications.



**StressMarq Biosciences**

June 15, 2016:

## Product Citations

---

FOR REFERENCE USE ONLY

93/100 Bioz Rating

11 Citations 2 Images

- Export this page of
- Export first 100 cita
- Export first 100 cita

**Bioz AI Summary:** The **HSP90 alpha ELISA Kit** from **StressMarq** has been extensively utilized across vario

Search term

Techniques

- Search for Techniques
- Enzyme-linked Immunosorbent Assay (ELISA)
- Protein Extraction (PROT-EX)
- Sandwich ELISA (ELISA-S)
- Spectrophotometry (SP-PH)

Journals

- Search for Journals
- American Journal of Physiology - Regulatory Integrative and Comparative Physiology
- Amyotrophic lateral sclerosis & frontotemporal degeneration
- Animal Production Science
- Europe PMC
- Gene

Dates

**Relationships between ambient temperature and diurnal variation in physiological responses to acute handling**  
 Pierre Danubian, Sarah E Polekoff, ..., Katherine L Buchanan  
 General and comparative endocrinology | 2025 Jun 10 | 40505874 | [Read Article](#) | [Bioz AI Brief](#)

**"Bioz AI Summary HSP90 alpha ELISA Kit from StressMarq was integral in measuring Hsp90 levels in house**  
 shock proteins. This helped elucid..." [More...](#)

[ELISA](#) [INTER-A](#)

Cited 1 time

**Prenatal Acoustic Signals Influence Nestling Heat Shock Protein Response to Heat and Heterophil-to-Lym**  
 Eve Udino, Anaïs Pessato, .... Mvlene M Mariette



FOR REFERENCE USE ONLY