



ELISA PRODUCT INFORMATION & MANUAL

HSP27 [p Ser15] *NBP2-62151*

Enzyme-linked Immunosorbent Assay for
quantitative detection of Human HSP27 [p Ser15].

For research use only.

Not for diagnostic or therapeutic procedures.

www.novusbio.com - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - technical@novusbio.com

Novus kits are guaranteed for 6 months from date of receipt

Product Manual

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Novus Biologicals is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Novus Biologicals' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Novus Biologicals, within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Novus Biologicals products and product applications are covered by US and foreign patents and patents pending. Novus Biologicals is a trademark of Novus Biologicals

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

Product Manual

TABLE OF CONTENTS



Please read
entire booklet
before
proceeding with
the assay.



Please contact
Novus
Biologicals
Technical
Support if
necessary.

Introduction	2
Principle	2
Materials Supplied.....	3
Storage	4
Other Materials Needed	4
Reagent Preparation	4
Sample Handling.....	6
Assay Procedure.....	8
Calculation of Results	9
Typical Results.....	10
Performance Characteristics	10
References.....	13

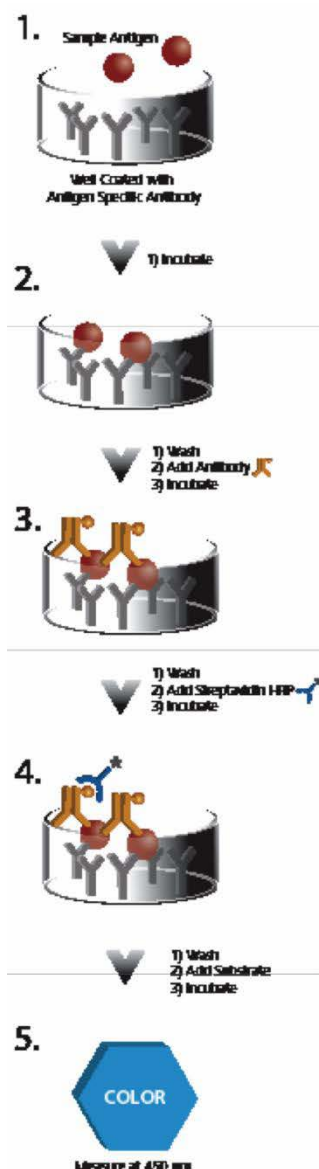
Product Manual

INTRODUCTION

Hsp27 is one of the most common members of the highly conserved and ubiquitously expressed family of small heat shock proteins (sHsp), which also includes α B-crystallin¹. It is characterized by a conserved C-terminal α -crystallin domain consisting of two anti-parallel β -sheets that promote oligomer formation required for its primary chaperone function as an inhibitor of irreversible protein aggregation². Hsp27 oligomerization is modulated by post-translational phosphorylation of Hsp27 at three serine residues, Ser15, Ser78, and Ser82, by a variety of protein kinases including MAPKAPK-3, PKA α , p70S6K, PKD I, and PKC δ ^{3,4}. Hsp27 has been shown to inhibit actin polymerization by binding of unphosphorylated Hsp27 monomers to actin intermediate filaments⁵. Anti-apoptotic functions of Hsp27 have also been identified through interactions with DAXX7, activation of Akt, and inhibition of apoptosome formation⁶⁻⁸. Evidence suggests altered expression of Hsp27 is implicated in the pathogenesis of breast, ovarian, and prostate cancer⁹.

PRINCIPLE

1. Samples and standards are added to wells coated with a monoclonal antibody specific for Hsp27. The plate is then incubated.
2. The plate is washed, leaving only bound Hsp27 on the plate. A yellow solution of antibody, specific for Hsp27 phosphorylated at Ser15, is then added. This binds the Hsp27 [p Ser15] captured on the plate. The plate is then incubated.
3. The plate is washed to remove excess antibody. A blue solution of HRP conjugate is added to each well, binding to the antibody, which is attached to the Hsp27 [p Ser15]. The plate is again incubated.
4. The plate is washed to remove excess HRP conjugate. TMB Substrate solution is added. The substrate generates a blue color when catalyzed by the HRP.
5. Stop solution is added to stop the substrate reaction. The resulting yellow color is read at 450 nm. The amount of signal is directly proportional to the level of Hsp27 [p Ser15] in the sample.



Product Manual

MATERIALS SUPPLIED



Do not mix components from different kit lots or use reagents beyond the expiration date of the kit.



Protect from light



Stop solution is caustic. Keep tightly capped.

- 1. Assay Buffer 27**
100 mL
Phosphate buffered saline containing BSA
- 2. Phospho Hsp27 Standard**
0.25 mL
One vial containing 10,000 pg/mL of recombinant human Hsp27 [p Ser15]
- 3. 5X Extraction Reagent**
10 mL
- 4. Hsp27 Clear Microtiter Plate**
One Plate of 96 Wells
A plate of break-apart strips coated with a mouse monoclonal antibody specific for Hsp27
- 5. Wash Buffer Concentrate**
100 mL
Tris buffered saline containing detergents
- 6. Hsp27 [p Ser15] ELISA Antibody**
10 mL
A yellow solution of rabbit polyclonal antibody specific for Hsp27, phosphorylated at Ser15
- 7. Hsp27 [p Ser15] ELISA Conjugate**
10 mL
A blue solution of mouse anti-rabbit IgG conjugated to horseradish peroxidase
- 8. TMB Substrate**
10 mL, Catalog No. 80-0350
A solution of 3,3',5,5' tetramethylbenzidine (TMB) and hydrogen peroxide
- 9. Stop Solution 2**
10 mL, Catalog No. 80-0377
A 1N solution of hydrochloric acid in water
- 10. Hsp27 [p Ser15] Assay Layout Sheet**
1 each, Catalog No. 30-0248
- 11. Plate Sealer**
3 each, Catalog No. 30-0012

Product Manual



Reagents require separate storage conditions.



Avoid repeated freeze/ thaw cycles.

STORAGE

All components of this kit, except the standard and antibody, are stable at 4°C until the kit's expiration date. The standard must be stored at or below -20°C upon receipt. Store the antibody at 4°C for up to 2 weeks; for long term storage beyond 2 weeks, store at -20°C. Avoid repeated freeze thaws.

OTHER MATERIALS NEEDED

1. Deionized or distilled water.
2. Phenylmethylsulfonyl fluoride (PMSF)
3. Protease inhibitor cocktail (PIC)
4. Phosphatase inhibitor cocktail (PhIC)
5. Precision pipets for volumes between 5 μ L and 1,000 μ L.
6. Repeater pipet for dispensing 100 μ L.
7. Disposable beakers for diluting buffer concentrates.
8. Graduated cylinders.
9. A microplate shaker.
10. Lint-free paper for blotting.
11. Microplate reader capable of reading at 450 nm.
12. Graph paper for plotting the standard curve.



Bring all reagents to room temperature for at least 30 minutes prior to opening.

REAGENT PREPARATION

1. Wash Buffer

Prepare the wash buffer by diluting 50 mL of the supplied Wash Buffer Concentrate with 950 mL of deionized water. This can be stored at room temperature until the kit's expiration, or for 3 months, whichever is earlier.

2. Extraction Reagent

Prepare the Extraction Reagent by diluting 10 mL of the supplied 5X Extraction Reagent with 40 mL of deionized water. This can be stored at 4°C until the kit expiration, or for 3 months, whichever is earlier.



If inhibitors other than those recommended are used, the end-user is responsible for assay validation.

Product Manual

3. PIC, PhIC, and PMSF Addition

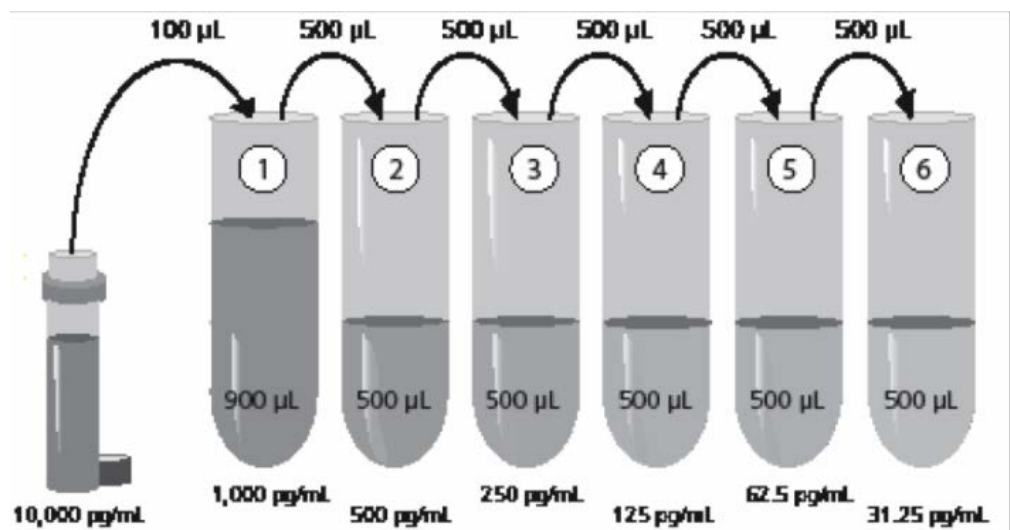
Immediately prior to use; PIC, PhIC, and PMSF must be added to the Extraction Reagent. Add 0.5 $\mu\text{L}/\text{mL}$ PIC, or equivalent concentration according to vendor's specification sheet. Add 10 $\mu\text{L}/\text{mL}$ PhIC, or equivalent concentration according to alternate vendor's specification sheet. Add PMSF to a final concentration of 1 mM.

Inhibitors must be freshly added to the Extraction Reagent to ensure optimal integrity of the samples. Each inhibitor treated buffer should incubate for 5-10 minutes at room temperature before it is used. Buffers treated with inhibitors should be used within 1 hour of preparation.

4. Preparation of phospho Hsp27 Standard Curve



Plastic tubes must be used for standard preparation.



Label six 12x75 mm polypropylene tubes #1 through #6. Pipet 900 μL of the assay buffer into tube #1. Pipet 500 μL of the assay buffer into tubes #2 through #6. Add 100 μL of the 10,000 pg/mL standard into tube #1. Vortex thoroughly. Add 500 μL of tube #1 to tube #2 and vortex thoroughly. Continue this for tubes #3 through #6.

Diluted standards should be used within 1 hour of preparation.

The concentrations of the standards are labeled above.

Product Manual

SAMPLE HANDLING



If buffers other than those provided are used, the end-user must determine the appropriate dilution and assay validation.



Samples should be stored at or below -20°C to avoid loss of bioactive analyte. Avoid repeated freeze/ thaw cycles.

This assay is suitable for measuring Hsp27 phosphorylated at Ser15 in cell lysates, tissue extracts, serum, and plasma. Prior to assay, frozen samples should be slowly brought to 4°C and centrifuged, if necessary, to isolate residual debris.

For cell lysates, a minimum 1:2 dilution of the 1X extraction reagent into the assay buffer is required to remove matrix interference of this buffer. Due to differences in sample types, number of cells, or total cellular protein concentration, samples may require greater dilution with the assay buffer to remove interference or to be read within the standard range. Users must determine the optimal sample dilutions for their particular experiments. Below are examples of sample recoveries with several different sample types. Note that % recovery was calculated based on linearity of samples.

Sample	Total cellular protein (mg/mL)	% Recovery	Recommended Dilution
HeLa Cells (human)	1.6	92%	1:800
Liver Microsomes (human)	1.0	103%	1:32
Serum (human)	N/A	99%	1:2
EDTA Plasma (human)	N/A	103-109%	1:64-1:512

Cell Lysate Preparation

1. Harvest cells and centrifuge at 250 x g (~1534 rpm) for 7 minutes at 4°C. Discard supernatant.
2. Resuspend pellet and wash with Hank's Balanced Salt Solution (without phenol red) or PBS.
3. Centrifuge at 250 x g (~1534 rpm) for 7 minutes at 4°C. Discard supernatant.
4. Resuspend pellet with Extraction Reagent plus inhibitors (see Reagent Preparation).
5. Vortex and incubate on ice for 30 minutes.
6. Centrifuge at 16,000 x g (~12,274 rpm) for 20 minutes at 4°C.
7. Place the supernatants into a clean tube.
8. The supernatants may be aliquoted and stored at or below -20°C or used immediately in the assay.



Add inhibitors to buffers prior to preparing samples

Product Manual

Serum and Plasma Preparation

1. Collect whole blood in either clotting tubes for serum or EDTA tubes for plasma.
2. Allow serum to clot for 30 minutes.
3. Centrifuge at 1000 x g for 15 minutes at 4°C.
4. Place supernatants into a clean tube.
5. The supernatant may be aliquoted and stored at or below -20°C, or used immediately in the assay.

Preparation of Tissue Extracts

1. Harvest tissue to be analyzed.
2. If necessary, tissues can be flash frozen and stored at -70°C. The extract can be prepared at a later time.
3. Calculate the amount of 1X Extraction Reagent that will be required. For each ~0.5cm³ piece of tissue, use 1mL of 1X Extraction Reagent.
4. Dilute an appropriate amount of 5X Extraction Reagent with cold (4°C) deionized or distilled water to generate the required volume of 1X Extraction Reagent. For example, if 5 mL of 1X Extraction Reagent were required, dilute 1 mL of the 5X Extraction Reagent with 4ml of cold deionized or distilled water.
5. Add protease and phosphatase inhibitors to the 1X Extraction Reagent.
6. Place the tissue in a mortar and add a sufficient volume of liquid nitrogen to cover the tissue.
7. Allow the liquid nitrogen to evaporate. The tissue should be thoroughly frozen.
8. Grind the frozen tissue to a powder with a pestle.
9. Add an appropriate volume of 1X Extraction Reagent supplemented with protease inhibitors to the processed tissue.
10. Continue to homogenize the tissue with the pestle until the tissue suspension is homogeneous.
11. Transfer the extract to a polypropylene tube and centrifuge at 21,000 x g for 10 minutes in a 4°C refrigerated microfuge.
12. Transfer the supernatant to a labeled polypropylene tube. The supernatant collected is the tissue extract, which is now ready for analysis using the Hsp27 [p Ser15] kit. The resulting pellet can be discarded.

Product Manual

13. Alternatively, the tissue extracts can be frozen at -70°C and assayed at a later date. It is recommended that a protein determination assay be performed and the extracts aliquoted into smaller volumes prior to storing at -70°C to avoid multiple freeze thaw cycles.



Bring all reagents to room temperature for at least 30 minutes prior to opening.



All standards, controls, and samples should be run in duplicate.



Pre-rinse each pipet tip with reagent. Use fresh pipet tips for each sample, standard, and reagent.



Pipet the reagents to the sides of the wells to avoid possible contamination



Prior to the addition of antibody, conjugate, and substrate, ensure there is no residual wash buffer in the wells. Remaining wash buffer may cause variation in assay results.

ASSAY PROCEDURE

Refer to the Assay Layout Sheet to determine the number of wells to be used.

Remove the wells not needed for the assay and return them, with the desiccant, to the mylar bag and seal. Store unused wells at 4°C .

1. Pipet 100 μL of the assay buffer into the S0 (0 pg/mL standard) wells.
2. Pipet 100 μL of Standards #1 through #6 to the bottoms of the appropriate wells.
3. Pipet 100 μL of the samples to the bottoms of the appropriate wells.
4. Seal the plate. Incubate for 1 hour shaking* at room temperature.
5. Empty the contents of the wells and wash by adding 400 μL of wash buffer to every well. Repeat 5 more times for a total of 6 washes. After the final wash, empty or aspirate the wells and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.
6. Pipet 100 μL of yellow antibody into each well except the blank.
7. Seal the plate. Incubate for 1 hour shaking* at room temperature.
8. Wash as above (Step 5).
9. Add 100 μL of blue conjugate to each well except the blank.
10. Seal the plate. Incubate for 30 minutes on a plate shaker* at room temperature.
11. Wash as above (Step 5).
12. Pipet 100 μL of substrate solution into each well.
13. Incubate for 30 minutes shaking* at room temperature.
14. Pipet 100 μL of stop solution into each well.
15. After blanking the plate reader against the substrate blank, read optical density at 450 nm. If plate reader is not capable of

Product Manual

adjusting for the blank, manually subtract the mean OD of the substrate blank from all readings.

**Shaking is preferably carried out on a suitable plate or orbital shaker set at a speed to ensure adequate mixing of the contents of the wells. The optimal speed for each shaker will vary and may range from 120-700 rpm.*

CALCULATION OF RESULTS

Several options are available for the calculation of the concentration of Hsp27 [p Ser15] in the samples. We recommend that the data be handled by an immunoassay software package utilizing a weighted 4-parameter logistic (4PL) curve fitting program.

1. Calculate the average Net Optical Density (OD) bound for each standard and sample by subtracting the average Blank OD from the average OD bound:

$$\text{Average Net OD} = \text{Average OD} - \text{Average Blank OD}$$



Make sure to multiply sample concentrations by the dilution factor used during sample preparation.

2. Plot Net OD versus Concentration of Hsp27 [p Ser15] for the standards. Sample concentrations may be calculated off of Net OD values using the desired curve fitting.

Make sure to multiply sample concentrations calculated off the curve by the dilution factor used during sample preparation to get starting sample concentration.

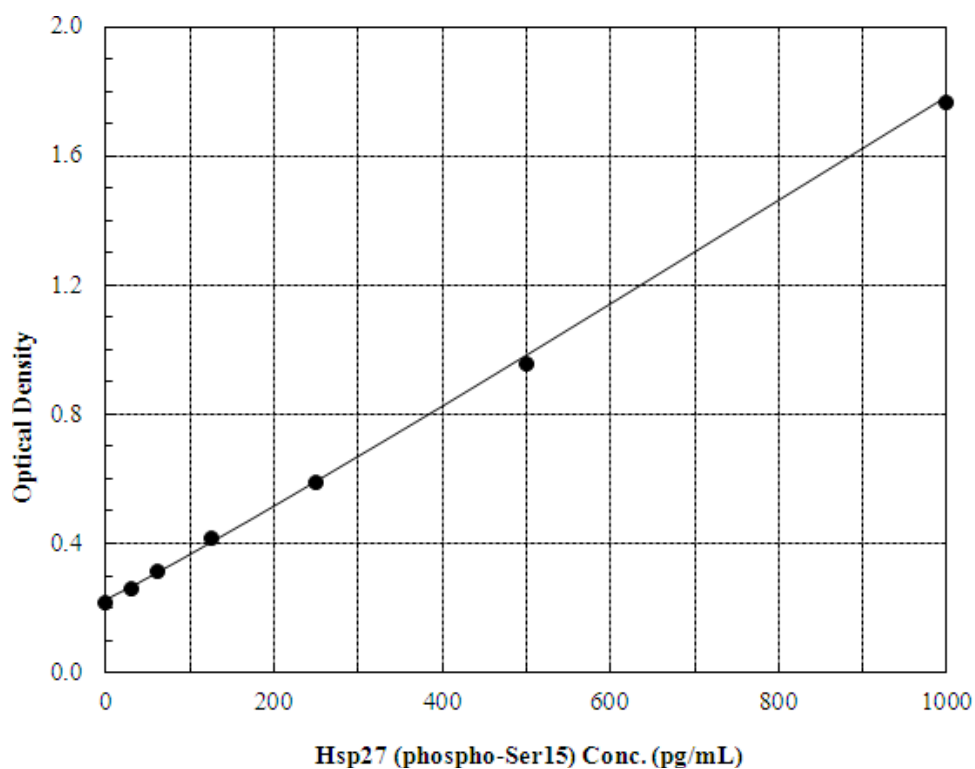
Samples with concentrations outside of the standard curve range will need to be re-analyzed using a different dilution.

Product Manual

TYPICAL RESULTS

The results shown below are for illustration only and should not be used to calculate results from another assay.

Sample	Net OD	Hsp27 (phospho-Ser15) (pg/mL)
S0	0.217	0
S1	1.768	1000
S2	0.955	500
S3	0.590	250
S4	0.417	125
S5	0.312	62.5
S6	0.262	31.25
Unknown 1	0.937	478
Unknown 2	0.309	62



PERFORMANCE CHARACTERISTICS

Specificity

The cross reactivities for a number of related compounds were determined by diluting cross reactants in the assay buffer at several concentrations. These samples were then measured in the assay.

Product Manual

Compound	Cross Reactivity
Hsp27 (non-phosphorylated)	0.885%
Hsp25	<0.016%
Hsp40	<0.016%
Hsp70	< 0.023%
Hsp90	< 0.016%
αB-Crystallin	< 0.016%
αA-Crystallin	< 0.016%

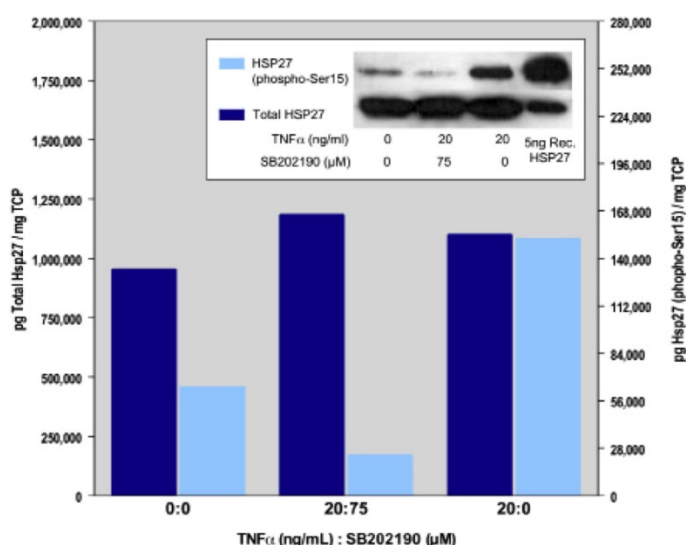
Hsp27 [p Ser15] has been shown to cross react at $\geq 100\%$. The percent cross reactivity calculation is relative to the phosphor Hsp27 standard used in the assay. The percent cross reactivity varies depending on the percent phosphorylation and protein concentration of the cross reactant used and has been shown to vary between lots.

Stimulation Experiments

This experiment was adapted from a protocol outlined in reference 10. Human HeLa cells were treated with 75 μM SB202190, a cell permeable inhibitor of p38 MAPK, for 1 hour at 37°C, followed by treatment with 20 ng/mL TNF- α , an inducer of Hsp27 phosphorylation, for 30 minutes at 37°C. Cells were washed 3 times in HBSS and

lysed in the 1X Extraction Reagent. Total cellular protein for each sample was determined using a BCA protein assay. Approximately 25 μg of total cellular protein was run on an 8-16% Tris-glycine gradient gel. Proteins were then transferred to a

nitrocellulose membrane and probed with antibodies against total Hsp27 and Hsp27 [p Ser15]. The same lysates were also diluted in the assay buffer and run in this kit, as well as a kit specific for total Hsp27, independent of phosphorylation.



Product Manual

Sensitivity

Sensitivity was calculated as the ratio of the mean OD plus 2 standard deviations of twenty-four replicates of the 0 pg/mL standard to the mean of twenty-four replicates of the lowest standard, multiplied by the concentration of that standard (31.25 pg/mL). This value was determined to be 10.15 pg/mL.

Linearity

A buffer sample containing Hsp27 [p Ser15] was serially diluted 1:2 in assay buffer and measured in the assay. The results are shown in the table below.

Dilution	Expected (pg/mL)	Observed (pg/mL)	Recovery (%)
Neat	---	918	---
1:2	459	449	98
1:4	230	195	85
1:8	115	93.6	81
1:16	57.4	57.1	100

Precision

Intra-assay precision was determined by assaying 20 replicates of three buffer controls containing Hsp27 [p Ser15] in a single assay.

pg/mL	%CV
449	4.0
110	9.3
48	14.2

Inter-assay precision was determined by measuring buffer controls of varying Hsp27 [p Ser15] concentrations in multiple assays over several days.

pg/mL	%CV
420	7.1
118	9.6
55	9.9

Product Manual

REFERENCES

1. Kappe, G., et al. (2003) Cell Stress Chaperones 8, 53-61.
2. Beck, F.X, et al. (2000) Am J Physiol Renal Physiol. 279, F203-F215.
3. Landry, J., et al. (1992) J Biol Chem. 267, 794-803.
4. Stokoe, D., et al. (1992) FEBS Lett. 313, 307-313.
5. Butt, E., et al. (2001) J Biol Chem. 276, 7108-7113.
6. Charette, S.J., et al. (2000) Mol Cell Biol. 20, 7602-7612.
7. Rane, M.J., et al. (2003) J Biol Chem. 278, 27828-27835.
8. Paul, C., et al. (2002) Mol Cell Biol. 22, 816-834.
9. Ciocca, D.R. and Calderwood, S.K. (2005) Cell Stress Chaperones 10, 86-103.
10. Park, K., et al. (2003) J Biol Chem. 37, 35272-35278.

Product Manual

Notes