



## **ELISA PRODUCT INFORMATION & MANUAL**

### **HSP27 [p Ser15] *NBP2-62152***

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

For research use only.

Not for diagnostic or therapeutic procedures.

**Human HSP27 [p Ser78] ELISA Kit**  
**Catalog # NBP2-62152**  
**96 Well Enzyme Immunoassay Kit**  
**For use with cell lysates, serum, and plasma**



All reagents, except standard, should be stored at 4°C. Store standard at -20°C.



Check our website for additional protocols, technical notes and FAQs.



For proper performance, use the insert provided with each individual kit received.

**Table of Contents**

<b><u>2</u></b>	<b><u>Introduction</u></b>
<b><u>2</u></b>	<b><u>Principle</u></b>
<b><u>3</u></b>	<b><u>Materials Supplied</u></b>
<b><u>4</u></b>	<b><u>Storage</u></b>
<b><u>4</u></b>	<b><u>Materials Needed but Not Supplied</u></b>
<b><u>5</u></b>	<b><u>Reagent Preparation</u></b>
<b><u>6</u></b>	<b><u>Sample Handling</u></b>
<b><u>7</u></b>	<b><u>Assay Procedure</u></b>
<b><u>8</u></b>	<b><u>Calculation of Results</u></b>
<b><u>9</u></b>	<b><u>Typical Results</u></b>
<b><u>10</u></b>	<b><u>Performance Characteristics</u></b>
<b><u>13</u></b>	<b><u>References</u></b>
<b><u>16</u></b>	<b><u>Limited Warranty</u></b>

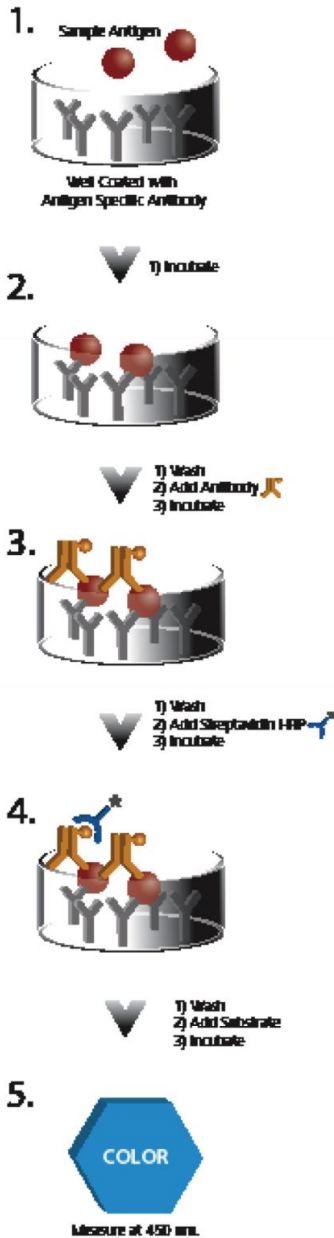
FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

## 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  $\alpha$ B-crystallin<sup>1</sup>. It is characterized by a conserved C-terminal  $\alpha$ -crystallin domain consisting of two anti-parallel  $\beta$ -sheets that promote oligomer formation required for its primary chaperone function as an inhibitor of irreversible protein aggregation<sup>2</sup>. 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 $\alpha$ , p70S6K, PKD I, and PKC $\delta$ <sup>3, 4</sup>. Hsp27 has been shown to inhibit actin polymerization by binding of unphosphorylated Hsp27 monomers to actin intermediate filaments<sup>5</sup>. Anti-apoptotic functions of Hsp27 have also been identified through interactions with DAXX7, activation of Akt, and inhibition of apoptosome formation<sup>6-8</sup>. Evidence suggests altered expression of Hsp27 is implicated in the pathogenesis of breast, ovarian, and prostate cancer<sup>9</sup>.

## 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 Ser78, is then added. This binds the Hsp27 (phospho-Ser78) 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 (phospho-Ser78). 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 (phospho-Ser78) in the sample.





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

## Materials Supplied

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 phospho Hsp27
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 (phospho-Ser78) EIA Antibody  
10 mL  
A yellow solution of mouse monoclonal antibody specific for Hsp27, phosphorylated at Ser78
7. Hsp27 (phospho-Ser78) EIA Conjugate  
10 mL  
A blue solution of mouse anti-rabbit IgG conjugated to horseradish peroxidase
8. TMB Substrate  
10 mL  
A solution of 3,3',5,5' tetramethylbenzidine (TMB) and hydrogen peroxide
9. Stop Solution 2  
10 mL  
A 1N solution of hydrochloric acid in water
10. Hsp27 (phospho-Ser78) Assay Layout Sheet  
1 each
11. Plate Sealer  
3 each



Protect substrate from prolonged exposure to light.



Stop solution is caustic. Keep tightly capped.



Reagents require separate storage conditions.

### Storage

All components of this kit, except the standard, are stable at 4°C until the kit's expiration date. The standard must be stored at or below -20°C.

### Materials Needed but Not Supplied

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  $\mu\text{L}$  and 1,000  $\mu\text{L}$ .
6. Repeater pipet for dispensing 100  $\mu\text{L}$ .
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.

## Reagent Preparation



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

### 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 45 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.

### 3. PIC, PhIC, and PMSF Addition

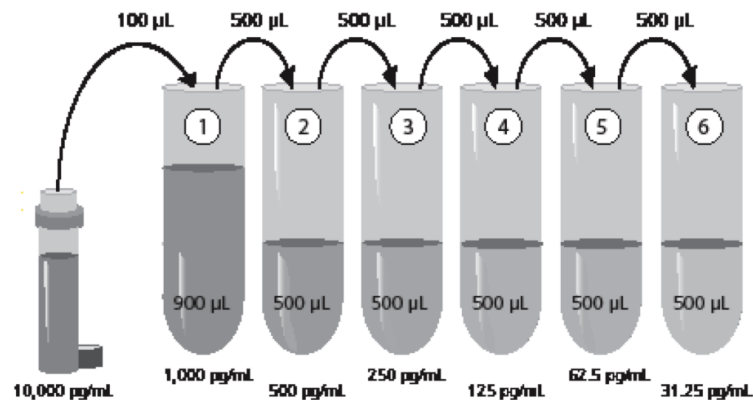
Immediately prior to use; PIC, PhIC, and PMSF must be added to the Extraction Reagent. Add 0.5 µL/mL PIC, or equivalent concentration according to vendor's specification sheet. Add 10 µL/mL PhIC, or equivalent concentration according to vendors specification sheet. Add PMSF to a final concentration of 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.



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



Samples must be stored frozen at or below -20°C to avoid loss of bioactive analyte. Repeated freeze/thaw cycles should be avoided.

## Sample Handling

This assay is suitable for measuring Hsp27 phosphorylated at Ser78 in cell lysates, 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)	1.55	89%	1:50
Serum (human)	N/A	99%	1:2
EDTA Plasma (human)	N/A	99%	1:100



Add inhibitors to buffers prior to preparing samples.

## 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.

## 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.

## Assay Procedure



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.

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  $\mu$ L of the assay buffer into the S0 (0 pg/mL standard) wells.
2. Pipet 100  $\mu$ L of Standards #1 through #6 to the bottoms of the appropriate wells.
3. Pipet 100  $\mu$ L 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  $\mu$ L 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  $\mu$ L 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  $\mu$ L of blue conjugate to each well except the blank.
10. Seal the plate. Incubate for 30 minutes on a plate shaker (~500 rpm) at room temperature.
11. Wash as above (Step 5).
12. Pipet 100  $\mu$ L of substrate solution into each well.
13. Incubate for 30 minutes shaking\* at room temperature.
14. Pipet 100  $\mu$ L 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 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 120-700rpm.



## Calculation of Results

Several options are available for the calculation of the concentration of Hsp27 (phospho-Ser78) in the samples. We recommend that the data be handled by an immunoassay software package utilizing a 4-parameter logistic curve fitting program. If data reduction software is not readily available, the concentrations can be calculated as follows:

1. Calculate the average Net OD for each standard and sample by subtracting the average blank OD from the average OD for each standard and sample.

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

2. Using linear graph paper, plot the average Net OD for each standard versus Hsp27 (phospho-Ser78) concentration in each standard. Approximate a straight line through the points. The concentration of the unknowns can be determined by interpolation.

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

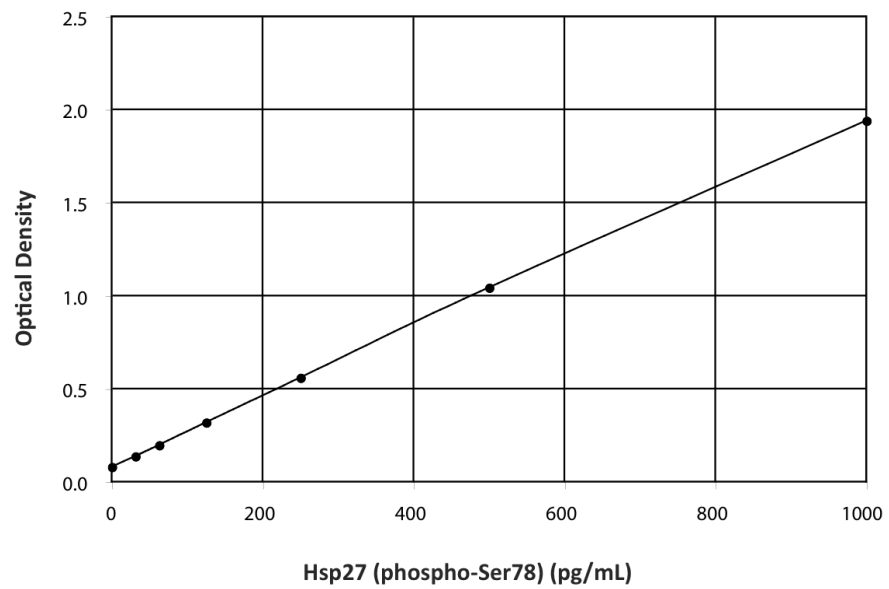


Make sure to multiply sample concentrations by the dilution factor used during sample preparation.

## 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-Ser78) (pg/mL)
S0	0.080	<b>0</b>
S1	1.940	<b>1000</b>
S2	1.044	<b>500</b>
S3	0.560	<b>250</b>
S4	0.319	<b>125</b>
S5	0.197	<b>62.5</b>
S6	0.137	<b>31.25</b>
Unknown 1	0.845	<b>396</b>
Unknown 2	0.185	<b>57</b>



## Performance Characteristics

### Specificity

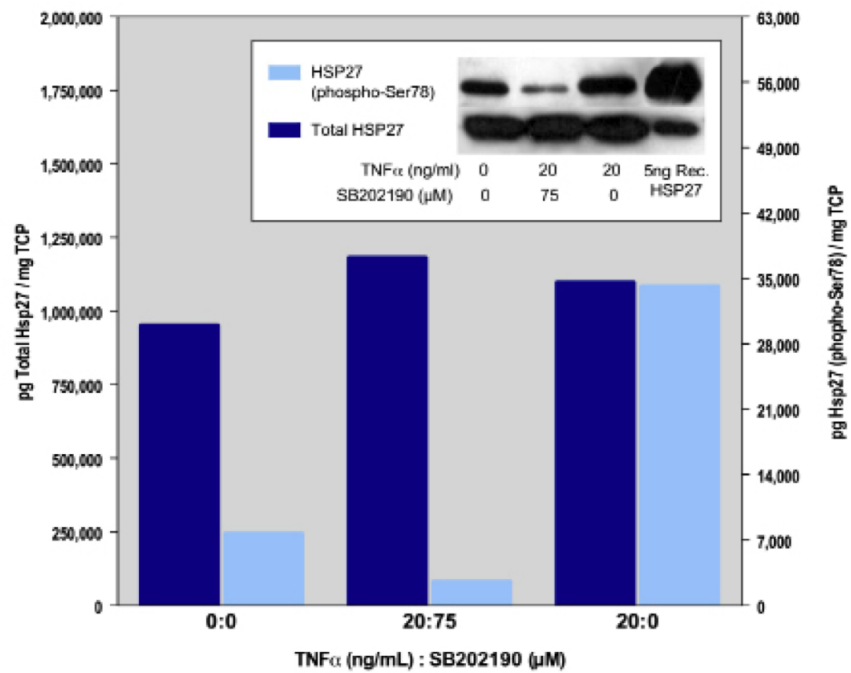
The cross reactivities for a number of related compounds were determined by diluting cross reactants in the assay buffer at several concentrations. These samples were then measured in the assay.

Compound	Cross Reactivity
Hsp27 (non-phosphorylated)	0.107%
Hsp25	<0.03%
Hsp40	<0.03%
Hsp70	< 0.03%
Hsp90	< 0.03%
$\alpha$ B-Crystallin	< 0.03%
$\alpha$ A-Crystallin	< 0.03%

Hsp27 (phospho-Ser78) has been shown to cross react at  $\geq 100\%$ . The percent cross reactivity calculation is relative to the phospho 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  $\mu$ M SB202190, a cell permeable inhibitor of p38 MAPK, for 1 hour at 37°C, followed by treatment with 20 ng/mL TNF- $\alpha$ , 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  $\mu$ g 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 (Cat. # SPA-803) and Hsp27 (phospho-Ser78) (Cat. # SPA-523PU). 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.



### 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 4.30 pg/mL.

### Linearity

A buffer sample containing Hsp27 (phospho-Ser78) 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	-----	783 pg/mL	-----
1:2	392 pg/mL	399 pg/mL	102 %
1:4	196 pg/mL	205 pg/mL	105 %
1:8	97.9 pg/mL	99.4 pg/mL	102 %
1:16	48.9 pg/mL	48.5 pg/mL	99 %

### Precision

**Intra-assay precision** was determined by assaying 20 replicates of three buffer controls containing Hsp27 (phospho-Ser78) in a single assay.

pg/mL	%CV
406	1.9
109	2.3
52	4.1

**Inter-assay precision** was determined by measuring buffer controls of varying Hsp27 (phospho-Ser78) concentrations in multiple assays over several days.

pg/mL	%CV
396	3.1
105	3.9
54	6.1

## 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.

## Notes

## Notes





MSDS (Material  
Safety Data Sheet)  
available online

### **Use of Product**

This product contains research chemicals. As such, they should be used and handled only by or under the supervision of technically qualified individuals. This product is not intended for diagnostic or human use.

### **Warranty**

Novus Biologicals makes no warranty of any kind, expressed or implied, which extends beyond the description of the product in this brochure, except that the material will meet our specifications at the time of delivery. Novus Biologicals makes no guarantee of results and assumes no liability for injuries, damages or penalties resulting from product use, since the conditions of handling and use are beyond our control.