



ELISA PRODUCT INFORMATION & MANUAL

GRP78/HSPA5 *NBP2-62145*

Enzyme-linked Immunosorbent Assay for quantitative detection of Human, Mouse, Rat GRP78/HSPA5.

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.



Carefully note the handling and storage conditions of each kit component.



Please contact Novus Biologicals Technical Support if necessary.

Background	2
Principle.....	3
Materials Supplied	4
Storage.....	5
Other Materials Needed.....	5
Sample Handling	6
Reagent Preparation.....	10
Assay Procedure	11
Calculation of Results	12
Typical Results	13
Performance Characteristics.....	14
References	16
Contact Information.....	18

Product Manual

BACKGROUND

The Novus Biologicals GRP78/HSPA5 Enzyme-Linked Immunosorbent Assay (ELISA) kit is a complete, competitive kit for the quantitative determination of GRP78/HSPA5 in cell lysates or serum. Please read the complete kit insert before performing this assay.

Glucose-regulated protein (GRP78/HSPA5) also known as binding immunoglobulin protein or BiP, is a resident molecular chaperone of the ER involved in the folding and assembly of proteins, transport of newly synthesized polypeptides across the ER membrane, regulation of calcium homeostasis and targeting misfolded proteins for degradation.¹⁻² GRP78 also regulates the transmembrane proteins, PERK, IRE1 and ATF6 by binding the N-terminal domains in the lumen of the ER preventing these signal transducers from initiating the unfolded protein response (UPR), an adaptive response to changes in the intracellular environment meant to restore normal ER function.³

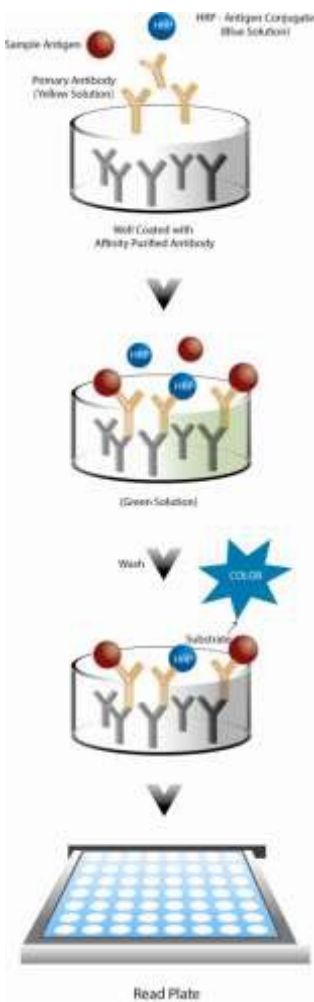
When protein misfolding occurs, exposed hydrophobic residues on the protein are bound by GRP78/HSPA5 preventing protein aggregation, further transit, and secretion. Intracellular stress such as glucose deprivation and viral infection can lead to a rapid accumulation of unfolded proteins.⁴ As unfolded proteins accumulate, more available GRP78/HSPA5 is required to bind the hydrophobic regions causing it to dissociate from the transmembrane signaling proteins thus initiating the UPR.

After dissociation from GRP78/HSPA5, activated PERK attenuates general translation to prevent further protein synthesis, ATF6 and IRE1 upregulate the production of ER folding and chaperone proteins including GRP78/HSPA5, and proteins that promote degradation of misfolded proteins through ER-associated protein degradation (ERAD). The overexpression of GRP78/HSPA5 as a result of UPR activation is believed to contribute to the pathology of metabolic disease, inflammatory disease, neurodegenerative disorders, and cancer.⁵⁻⁸

Product Manual

PRINCIPLE

1. Standards and samples are added to wells coated with a donkey anti-sheep IgG antibody. A solution of goat polyclonal antibody to GRP78/HSPA5 is also added to each well and the plate is incubated.
2. A solution of GRP78/HSPA5 conjugated to horseradish peroxidase is then added to each well and incubated. During this incubation the antibody binds, in a competitive manner, the GRP78/HSPA5 in the sample or conjugate. The plate is washed leaving only bound GRP78/HSPA5.
3. TMB substrate solution is added. An HRP-catalyzed reaction generates a blue color in the solution.
4. Stop solution is added to stop the substrate reaction. The resulting yellow color is read at 450 nm. The amount of signal is inversely proportional to the level of GRP78/HSPA5 in the sample.



Product Manual

MATERIALS SUPPLIED

1. Donkey anti-Sheep IgG Microtiter Plate

One plate of 96 wells

A clear plate of break-apart strips coated with a Donkey anti-sheep polyclonal antibody

2. Assay Buffer 28

50 mL

Tris buffered saline containing BSA and detergents

3. GRP78/HSPA5 Conjugate

5 mL

A blue solution of GRP78/HSPA5 conjugated to horseradish peroxidase

4. GRP78/HSPA5 Antibody

5 mL

A yellow solution of goat polyclonal antibody specific for GRP78/HSPA5

5. GRP78/HSPA5 Standard

Two vials each containing 4.5 µg of recombinant GRP78/HSPA5

6. Extraction Reagent #2

10 mL

5x Tris buffered solution containing detergent

7. Wash Buffer Concentrate

100 mL

20x Tris buffered saline containing detergents

8. TMB Substrate

2x10 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. GRP78/HSPA5 Assay Layout Sheet

1 each

11. Plate Sealer

3 each



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



Protect substrate from prolonged exposure to light.



Stop solution is caustic. Keep tightly capped.

Product Manual

STORAGE

All kit components should be stored at 4°C except the GRP78/HSPA5 Conjugate which needs to be stored at -80°C upon receipt. Avoid repeated freeze thaw cycles of the GRP78/HSPA5 Conjugate. Shipping conditions may not reflect storage conditions.

OTHER MATERIALS NEEDED

1. Deionized or distilled water
2. Phenylmethylsulfonyl fluoride (PMSF)
3. Protease Inhibitor Cocktail (PIC)
4. Precision pipets for volumes between 5 μ L and 1,000 μ L
5. Repeater pipet for dispensing 100 μ L
6. Disposable beakers for diluting buffer concentrates
7. Graduated cylinders
8. A microplate shaker
9. Adsorbent paper for blotting
10. Microplate reader capable of reading a 450 nm
11. Software for extrapolating sample values from optical density readings utilizing a four parameter logistic curve fit.



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



Sample handling procedures should be completed prior to reagent preparation.

SAMPLE HANDLING

The assay is suitable for the measurement of GRP78/HSPA5 in cell lysates and human serum. Prior to assay, frozen samples should be brought to 4°C and centrifuged, if necessary, to remove residual debris.

A minimum dilution of 1:5 in assay buffer is required to remove matrix and or lysis buffer interference with most cell lysate samples. It is necessary to validate the use of any lysis buffer other than that provided in this kit.

Due to variation in samples, a different dilution may be required. Users must determine the optimal dilutions for their particular experiments.

Linearity

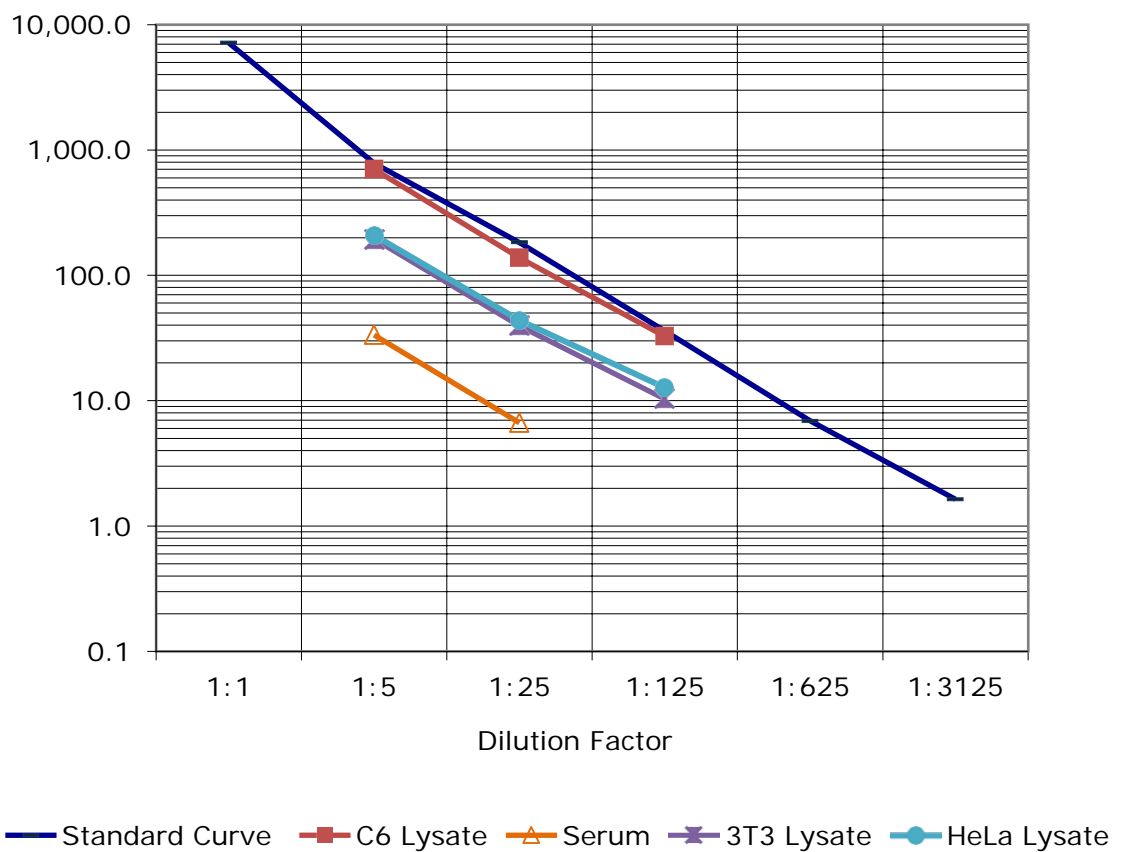
The minimum required dilution for common lysates and serum was determined by serially diluting samples into the provided assay buffer and identifying the dilution at which linearity was observed.

% Dilutional Linearity				
Dilution	HeLa Lysate	C6 Lysate	3T3 Lysate	Serum
Neat	---	---	---	---
1:5	96	101	99	101
1:25	100	100	100	100
1:125	116	118	103	---

Product Manual

Parallelism

To assess parallelism, HeLa, 3T3, and C6 lysates and serum were serially diluted into assay buffer and run in the assay. The GRP78/HSPA5 concentration in each sample was assigned using the standard curve. Assigned concentrations were plotted as a function of sample dilution. Parallelism of the curves demonstrates that the antigen binding characteristics are similar enough to allow the accurate determination of native analyte levels in diluted samples of human, mouse, and rat origin.



Product Manual

Spike and Recovery

After diluting each individual sample to read within the dynamic range of the assay, recombinant GRP78/HSPA5 was spiked at high, medium and low concentrations. Endogenous GRP78/HSPA5 was subtracted from the spiked values and the recovery in each of the spiked matrices was compared to the recovery of identical spikes in the assay buffer. The percent recovery at the three concentrations is indicated below for each matrix.

Sample Matrix	Dilution	Spike Concentration	Recovery of Spike
Hela Cell Lysate	1:25	1000 ng/mL	109%
		300 ng/mL	108%
		13 ng/mL	90%
C6 Cell Lysate	1:125	1000 ng/mL	82%
		300 ng/mL	93%
		13 ng/mL	101%
3T3 Cell Lysate	1:25	1000 ng/mL	94%
		300 ng/mL	97%
		13 ng/mL	98%
Serum	1:5	1000 ng/mL	120%
		300 ng/mL	110%
		13 ng/mL	113%

Product Manual

Cell Lysate Preparation

1. Centrifuge at 1700 x g for 10 minutes at room temperature to pellet cells and/or cellular debris.
2. Prepare Extraction Reagent
 - a. Calculate the amount of 1X Extraction Reagent that will be required. For every 1×10^6 to 1×10^7 cells, use 1 mL of 1X Extraction Reagent.
 - b. Dilute an appropriate amount of 5X Extraction Reagent with cold deionized or distilled water to generate the required volume of 1X Extraction Reagent.
 - c. Add 0.5 μ L of protease inhibitor cocktail (PIC) per mL of 1X Extraction Reagent and add PMSF to a final concentration of 1mM. Inhibitors must be added fresh just prior to lysis. Extraction Reagent containing inhibitors should not be stored for later use.
3. Resuspend cell pellet in 1X Extraction Reagent with inhibitors and incubate on ice for 30 minutes. Vortex occasionally.
4. Pellet cellular debris via centrifugation at 10,000 x g for 10 minutes in a 4°C refrigerated centrifuge.
5. Collect the supernatant (lysate) without disturbing the cell pellet. Divide the lysates into aliquots and store at or below -20°C, or use immediately in the assay.
6. Avoid repeated freeze-thaw cycles.

Serum Preparation

1. Collect whole blood in appropriate serum tubes.
2. Blood samples should be incubated in an upright position at room temperature for 30-45 min (no longer than 60 min) to allow clotting.
3. Centrifuge at 1000 x g for 15 minutes at room temperature.
4. Remove serum to a clean plastic tube.
5. Divide serum into aliquots and store at or below -20°C, or use immediately in the assay.
6. Avoid repeated freeze-thaw cycles.

REAGENT PREPARATION

1. Wash Buffer

Prepare Wash buffer by diluting 30 mL of the supplied Wash Buffer concentrate with 570 mL of deionized water. Store the diluted wash buffer at room temperature. Diluted wash buffer should be used within 3 months.

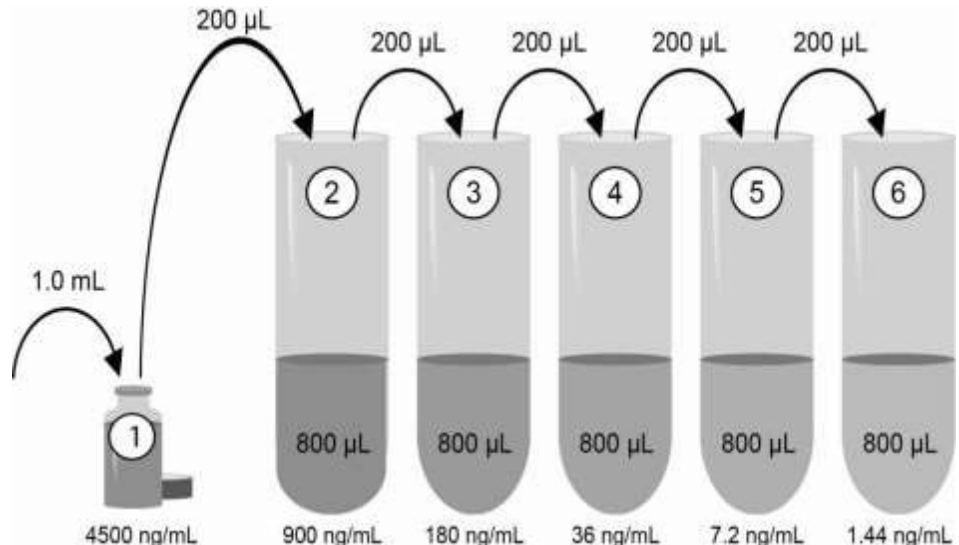
2. GRP78/HSPA5 Standard

Allow the GRP78/HSPA5 standard to warm to room temperature. Reconstitute one vial of GRP78/HSPA5 standard with 1.0 mL Assay Buffer 28 for a 4500ng/mL standard concentration. Vortex thoroughly, wait 5 minutes and vortex again prior to use. Label this vial #1.

Label five 12 x 75mm tubes #2 through #6. Pipet 800 μ L of assay buffer into each tube. Remove 200 μ L from the reconstituted standard vial and add to tube #2 and vortex thoroughly. Remove 200 μ L from tube #2 and add to tube #3. Vortex thoroughly. Continue this for tubes #4 through #6.



Polypropylene tubes may be used for standard preparation. Avoid polystyrene.



Diluted standards should be used immediately. Do not store remaining standard.

All other kit components should be brought to room temperature prior to use in the assay.

ASSAY PROCEDURE



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



All standards and samples should be run in duplicate.



Do Not Wash wells after step 5.



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

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 μ L of Assay Buffer 28 into the B₀ (0 ng/mL standard) wells.
2. Pipet 100 μ L of standards #1 through #6 to the bottom of the appropriate wells.
3. Pipet 100 μ L of the samples into the appropriate wells.
4. Pipet 50 μ L of yellow Antibody solution into each well, except the blanks.
5. Seal the plate. Incubate for 1 hour with mixing on a plate shaker at room temperature*. **Do not wash at this step.**
6. Add 50 μ L of blue conjugate to each well except the blank. Seal the plate. Incubate for 1 hour with mixing on a plate shaker at room temperature.
7. Empty the contents of the wells and wash with an automated washer by adding 300 μ L of 1X Wash Buffer to every well. Repeat 3 more times for a total of 4 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.
8. Pipet 200 μ L of TMB solution into each well.
9. Seal the plate. Incubate for 30 minutes with shaking on a plate shaker at room temperature.
10. Pipet 50 μ L of Stop Solution into each well.
11. After blanking the plate reader against the substrate, read optical density at 450 nm. If the plate reader is not capable of adjusting for the blank, manually subtract the mean OD of the substrate blank from all readings.

* **Note:** The optimal speed for each shaker will vary and may range from 120-700 rpm. The speed must be set to ensure adequate mixing of the wells, but not so vigorously that the contents of the wells splash out and contaminate other wells.



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

CALCULATION OF RESULTS

Several options are available for the calculation of the concentration of GRP78/HSPA5 in the samples. We recommend that the data be handled by an immunoassay software package utilizing a 4 parameter logistic (4PL) curve fitting program. The concentration of GRP78/HSPA5 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. Calculate the binding of each pair of standard wells as a percentage of the maximum binding wells (B_0), using the following formula:

$$\text{Percent Bound} = \frac{\text{Net OD}}{\text{Net } B_0 \text{ OD}} \times 100$$

3. Plot the Percent Bound (B/B_0) versus concentration of GRP78/HSPA5 for the standards. Approximate a straight line through the points. The concentration of GRP78/HSPA5 in the unknowns can be determined by interpolation.

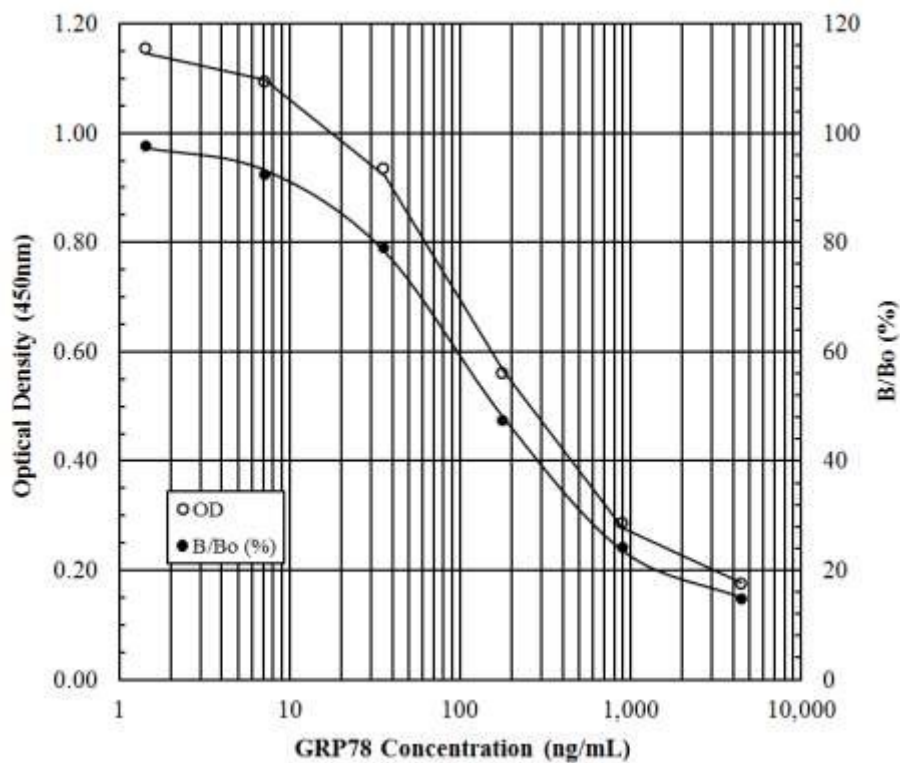
Samples with concentrations outside of the standard curve range will need to be re-analyzed using a different dilution.

Product Manual

TYPICAL RESULTS

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

Sample	Average Net OD	Percent Bound	GRP78/HSPA5 (ng/mL)
Blank	0.044	---	---
B ₀	1.182	100%	0
S1	0.174	14.7%	4500
S2	0.285	24.1%	900
S3	0.560	47.4%	180
S4	0.933	79.0%	36
S5	1.092	92.4%	7.2
S6	1.120	97.7%	1.44



Product Manual

PERFORMANCE CHARACTERISTICS

Specificity

The cross reactivities for related compounds were determined by diluting the cross reactant in the kit assay buffer at a concentration of ten times the high standard then measuring in the assay.

Analyte	Cross Reactivity
GRP94	≤0.01%
PDI	≤ 0.01%
Calreticulin	≤ 0.01%

Sensitivity

The sensitivity or limit of detection of the assay is 8.4 ng/mL, determined by interpolation at 2 standard deviations away from the mean signal of 24 replicates of zero standard using data from 8 standard curves.

Interference

Detergents commonly used in lysis buffers were analyzed for interference in the assay and the tolerance was determined.

Detergent	Assay Tolerance	
	(%)	molarity
SDS	≤ 0.03	≤ 1 mM
Tx100	≤ 1.0	≤ 15 mM
NP-40	≤ 6.0	≤ 98 mM
Sodium Deoxycholate	≤ 0.06	≤ 1 mM
Zwittergent	≤ 0.1	≤ 3.4 mM

Product Manual

Intra-assay precision was determined by assaying 24 replicates of two buffer controls containing GRP78/HSPA5 in a single assay.

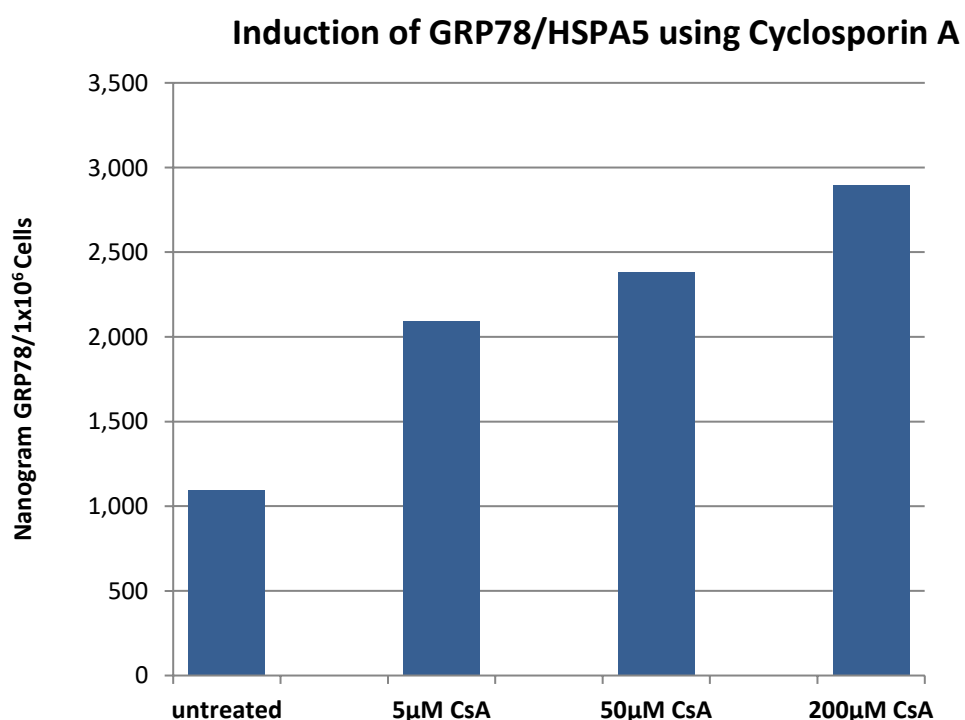
ng/mL	%CV
373.2	4.5
12.5	14.1

Inter-assay precision was determined by measuring buffer controls of varying GRP78/HSPA5 concentrations in multiple assays over several days.

ng/mL	%CV
352.8	4.8
13.0	13.1

Experimental Evaluation

HeLa cells were treated for 5 hours with 5 μ M, 50 μ M or 200 μ M concentrations of cyclosporine A. Following the incubation, treated and untreated cells were lysed with Extraction Reagent #2 as described above. The lysates were then evaluated in the assay. This data agrees with the results reported in the literature, low concentrations of cyclosporine A lead to the induction of GRP78/HSPA5 in HeLa cells (Paslaru, L. et al. (1994)) and supports the claim that the Novus ELISA is able to detect and quantitate changes in native levels of GRP78/HSPA5.



Product Manual

REFERENCES

1. Liu, C. (2003). The unfolded protein response. *Journal of Cell Science*, 116:1861-18662.
2. Schroder, M. (2006). The unfolded protein response. *Molecular Biotechnology*, 34:279-290.
3. Herbert, D et.al. (2007). In and out of the ER: protein folding, quality control, degradation, and related human diseases. *Physiological Reviews*, 87:1377-1408.
4. Xu, C. et.al. (2005). Endoplasmic reticulum stress: cell life and death decisions. *The Journal of Clinical Investigation*, 115(10):2656-2664.
5. Wang, S. et.al. (2012). The impact of the unfolded protein response on human disease. *The Journal of Cell Biology*, 197(7):587-867.
6. Wang, M. et.al. (2009). Role of the unfolded protein response regulator GRP78/BiP in development, cancer, and neurological disorders. *Antioxidants and Redox Signaling*, 11(9):2307-2316.
7. Lee, A. (2007). GRP78 induction in cancer: therapeutic and prognostic implications. *Cancer Research*, 67:3496-3499.
8. Zheng, H. et.al. (2008). Overexpression of GRP78 and GRP94 are markers for aggressive behavior and poor prognosis in gastric carcinomas. *Human Pathology*, 39:1042-1049.
9. Paslaru, L. et al. (1994). GRP78 induction by cyclosporine A in human HeLa cells. *FEBS*, 350:304-308.

Product Manual

NOTES

Product Manual

NOTES