



IMMUNOSET® HSP25 (rodent) ELISA development set

Catalog number: ADI-960-075
Reagents for 5 x 96 wells

This IMMUNOSET® HSP25 (rodent) ELISA development set contains the basic components for the development of an Hsp25 immunometric enzyme-linked immunosorbent assay (ELISA). Each kit contains sufficient reagents for five 96-well plates.

This kit has been validated for use with cell lysates. Additional sample types will require validation by the user. Visit www.enzolifesciences.com for tips and frequently asked questions.

Introduction

Rodent Hsp25, human Hsp27, and α B-crystallin are part of a diverse family of small heat shock proteins (sHsps) that range in size from 14-45 kDa¹. sHsps are constitutively expressed, and upregulated in response to stress such as heat shock. They function as chaperone-like proteins by binding unfolded polypeptides and preventing uncontrolled protein aggregation. Data indicates that Hsp25 is a dynamic tetramer of tetramers with a unique ability to refold and reassemble into its active quaternary structure after denaturation². Hsp25 helps facilitate the glutathione-redox cycle by enhancing glutathione utilization and maintaining the cellular glutathione pool in favor of the reduced states³.

References:

1. Narberhaus, F. (2002) *Microbiol Mol Biol Rev.* **66**, 64-93.
2. Ehrnsperger, M., *et al.* (1999) *J Biol Chem.* **274**, 14867-14874.
3. Baek, S.H., *et al.* (2000) *J Cell Physiol* **183**, 100-107.

Materials Provided

1. Hsp25 Capture Antibody
One vial containing lyophilized Hsp25 monoclonal antibody, Component #80-1942
2. Hsp25 Standard
One vial containing 312.5 ng lyophilized recombinant Hsp25 protein, Component #80-1943
3. Hsp25 Detection Antibody
One vial containing lyophilized Hsp25 biotinylated polyclonal antibody, Component #80-1944
4. SA-HRP
One vial containing 12.5 μ g lyophilized streptavidin conjugated to horseradish peroxidase, Component #80-1896

Materials Needed but not Supplied

1. RIPA Cell Lysis Buffer, Cat. #80-1284, or similar
2. 96-well high-binding polystyrene microtiter plates, Cat. #80-1930, or similar
3. Precision pipets
4. Microplate reader capable of reading at 450 nm
5. Microplate shaker
6. Phosphate buffered saline (PBS)[†]
7. Tween[®]-20^{*†}
8. Bovine Serum Albumin (BSA)[†]
9. 3,3',5,5' tetramethylbenzidine (TMB) solution, Cat. #80-1805 or similar[†]
10. 1N hydrochloric acid, such as Stop Solution 2, Cat. #80-1804[†]
11. Sucrose
[†]IMMUNOSET® Buffer Pack, Cat. #ADI-950-003
^{*}Tween is a registered trademark of ICL Americas

Buffer Formulations

1. Coating Buffer
10 mM sodium phosphate, 15 mM NaCl, pH 7.4
2. Blocking Buffer
10 mM sodium phosphate, 15 mM NaCl, 1.0% BSA, 1% sucrose, pH 7.4
3. Assay Buffer
100 mM sodium phosphate, 150 mM NaCl, 1.0% BSA, 0.1% Tween-20, pH 7.4
4. Wash Buffer
10 mM sodium phosphate, 15 mM NaCl, 0.1% Tween-20, pH 7.4

Plate Coating

1. Reconstitute Hsp25 Capture Antibody with 250 μ L deionized water for a 250x stock. Use immediately, or make aliquots and freeze at -20°C for up to 3 months. For prolonged storage, aliquot and freeze at -80°C. Avoid repeated freeze/thaw cycles.
2. Dilute the stock 1:250 in Coating Buffer. Immediately dispense into 96-well microtiter plates using 100 μ L of the diluted capture antibody per well. Seal the plate and incubate overnight at room temperature.
3. Aspirate each well to remove coating solution. Immediately add 200 μ L Blocking Buffer per well. Seal the plate and incubate for at least 1 hour.
4. Aspirate each well to remove blocking solution. Plates may be used immediately or dried and stored with desiccant at 4°C.

Reagent Preparation

1. Recombinant Hsp25 Standard
Reconstitute vial contents with 250 μ L deionized water for a 1250 ng/mL (50x) stock. Aliquot and store at -20°C for up to 3 months. For prolonged storage, aliquot and freeze at -80°C. Avoid repeated freeze/thaw cycles.

The recommended standard curve range is 25 ng/mL to 0.78 ng/mL, using 2-fold serial dilutions in Assay Buffer. Do not store diluted standard.
2. Hsp25 Detection Antibody
Reconstitute vial contents with 250 μ L deionized water for a 250x stock. Aliquot and store at -20°C for up to 3 months. For prolonged storage, aliquot and freeze at -80°C. Avoid repeated freeze/thaw cycles.

Dilute the stock 1:250 in Assay Buffer for a working solution. Do not store diluted antibody.
3. SA-HRP
Reconstitute vial contents with 250 μ L deionized water for a 1500x stock. Store at 4°C for up to 3 months. For prolonged storage, aliquot and freeze at -20°C. Avoid repeated freeze/thaw cycles.

Dilute the stock 1:1500 in Assay Buffer for a working solution. Do not store diluted conjugate.

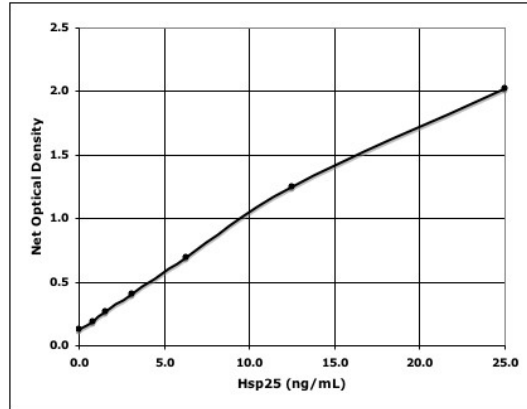
Assay Procedure

1. Pipet 100 μ L of Assay Buffer into the control (0 ng/mL standard) wells.
2. Pipet 100 μ L of standards and samples, prepared in Assay Buffer, to the bottom of the appropriate wells.
3. Seal the plate. Incubate for 1 hour on a plate shaker at room temperature.
4. Empty the contents of the wells and wash by adding 400 μ L of 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.
5. Pipet 100 μ L of the diluted detection antibody into each well, except the blank.
6. Seal the plate. Incubate for 1 hour on a plate shaker at room temperature.
7. Wash as above (Step 4).
8. Add 100 μ L of the diluted conjugate to each well except the blank.
9. Seal the plate. Incubate for 30 minutes on a plate shaker at room temperature.
10. Wash as above (Step 4).
11. Pipet 100 μ L of TMB solution into each well.
12. Seal the plate. Incubate for 30 minutes on a plate shaker at room temperature.
13. Pipet 100 μ L 1N HCl into each well.
14. 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.

Assay Performance

Typical Data

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



Sensitivity

The sensitivity, or limit of detection, of this assay is 0.38 ng/mL. It was determined by interpolation at 2 standard deviations above the mean signal at background, using data from 5 standard curves.

Specificity

This assay detects Hsp25 in cell lysates of rat and mouse origin. No cross reactivity is observed with human Cpn10, Hsp27, and Hsp40, rat Hsp32, or bovine α B Crystallin.

Dilutional Linearity

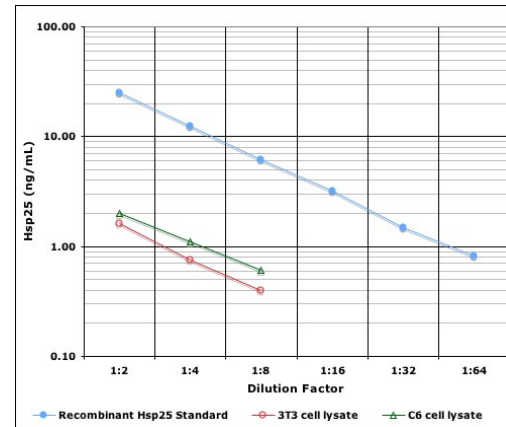
To determine possible interference from the sample matrix, the indicated sample types were serially diluted into assay buffer. The concentrations of Hsp25 were measured in the assay, and the results were analyzed to determine the range over which a linear response was obtained. These data may be used as a guideline to determine minimal recommended dilution (MRD) for similar samples.

Dilution Factor	C6 CL	3T3 CL
1:2	29%	30%
1:4	61%	78%
1:8	82%	102%
1:16	90%	94%
1:32	100%	100%

CL: Cell Lysate

Parallelism

Dose-response curves from cell lysates diluted into assay buffer (using the MRD) were compared to the recombinant Hsp25 standard curve. Parallelism indicates antibody-binding characteristics of the native and standard proteins are similar, allowing accurate determination of analyte.



Calculation of Results

Several options are available for the calculation of the relative levels of Hsp25 in samples. We recommend that the data be handled by an immunoassay software package utilizing a 4 parameter logistic curve-fitting program. For accuracy, please ensure that sample values fall within the standard range.

Accessory Reagent List		
Reagent	Quantity	Cat. #
IMMUNOSET [®] Buffer Pack	1 each of the following products: 80-1927, 80-1928, 80-1929, 80-1805, 80-1804	ADI-950-003
IMMUNOSET [®] Plate Pack	5 96-well clear microtiter plates & 5 plate sealers	80-1930
PBS Concentrate	120 mL	80-1927
BSA Solution (10%)	50 mL	80-1928
Tween-20 Solution (10%)	30 mL	80-1929
RIPA Cell Lysis Buffer 2	100 mL	80-1284
Wash Buffer Concentrate	100 mL	80-1287
SA-HRP	12.5 μ g/vial	80-1896

Storage

Store all components at 4°C. See page 3 for storage of reconstituted material.

Tips & Troubleshooting

- ✓ If buffers other than those recommended are used in the assay, the end-user must determine the appropriate dilution and assay validation.
- ✓ Pipet the reagents to the sides of the wells to avoid possible contamination.
- ✓ Pre-rinse each pipet tip with reagent. Use fresh pipet tips for each sample, standard, and reagent.
- ✓ Insufficient washing or residual wash buffer in the wells may cause variation in assay results.
- ✓ Bring all reagents to room temperature for at least 30 minutes prior to opening.
- ✓ All standards, controls, and samples should be assayed in duplicate.

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 use only. Not for use in diagnostic procedures. 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 Enzo Life Sciences, Inc. 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

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. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., 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.

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