

Version 2b Last updated 28 February 2023

ab233620

Mouse CD93

SimpleStep ELISA[®] Kit

For the quantitative measurement of CD93 in mouse serum, plasma, tissue supernatant, and tissue homogenate extract samples.

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Overview	1
2. Protocol Summary	2
3. Precautions	3
4. Storage and Stability	3
5. Limitations	4
6. Materials Supplied	4
7. Materials Required, Not Supplied	5
8. Technical Hints	5
9. Reagent Preparation	7
10. Standard Preparation	8
11. Sample Preparation	9
12. Plate Preparation	11
13. Assay Procedure	12
14. Calculations	14
15. Typical Data	15
16. Calibration	17
17. Typical Sample Values	17
18. Assay Specificity	24
19. Species Reactivity	24
20. Troubleshooting	25
Technical Support	26

1. Overview

CD93 *in vitro* SimpleStep ELISA® (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of CD93 protein in mouse serum, plasma, tissue supernatant, and tissue homogenate extract samples

The SimpleStep ELISA® employs an affinity tag labeled capture antibody and a reporter conjugated detector antibody which immunocapture the sample analyte in solution. This entire complex (capture antibody/analyte/detector antibody) is in turn immobilized via immunoaffinity of an anti-tag antibody coating the well. To perform the assay, samples or standards are added to the wells, followed by the antibody mix. After incubation, the wells are washed to remove unbound material. TMB Development Solution is added and during incubation is catalyzed by HRP, generating blue coloration. This reaction is then stopped by addition of Stop Solution completing any color change from blue to yellow. Signal is generated proportionally to the amount of bound analyte and the intensity is measured at 450 nm. Optionally, instead of the endpoint reading, development of TMB can be recorded kinetically at 600 nm.

CD93 is a transmembrane glycoprotein receptor for C1q, MBL2, and SPA. It is involved in the immune response and in hematopoietic and vascular development. Elevated CD93 levels are seen during peritonitis, sepsis, and rheumatoid arthritis. Altered CD93 serum levels correlates with a decreased risk of coronary artery disease and myocardial infarction.

2. Protocol Summary

Prepare all reagents, samples, and standards as instructed



Add 50 μ L standard or sample to appropriate wells



Add 50 μ L Antibody Cocktail to all wells



Incubate at room temperature for 1 hour



Aspirate and wash each well three times with 350 μ L 1X Wash Buffer
PT



Add 100 μ L TMB Development Solution to each well and incubate
for 10 minutes.



Add 100 μ L Stop Solution and read OD at 450 nm

3. Precautions

Please read these instructions carefully prior to beginning the assay.

- All kit components have been formulated and quality control tested to function successfully as a kit.
- We understand that, occasionally, experimental protocols might need to be modified to meet unique experimental circumstances. However, we cannot guarantee the performance of the product outside the conditions detailed in this protocol booklet.
- Reagents should be treated as possible mutagens and should be handled with care and disposed of properly. Please review the Safety Datasheet (SDS) provided with the product for information on the specific components.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas.
- All biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

4. Storage and Stability

Store kit at +4°C immediately upon receipt. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.

Refer to list of materials supplied for storage conditions of individual components.

5. Limitations

- Assay kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted.

6. Materials Supplied

Item	Quantity	Storage Condition
Mouse CD93 Capture Antibody 10X	600 µL	+4°C
Mouse CD93 Detector Antibody 10X	600 µL	+4°C
Mouse CD93 Lyophilized Recombinant Protein	2 Vials	+4°C
Antibody Diluent 4BR	6 mL	+4°C
Wash Buffer PT 10X	20 mL	+4°C
Cell Extraction Buffer PTR 5X	10 mL	+4°C
Cell Extraction Enhancer Solution 50X	1 mL	+4°C
TMB Development Solution	12 mL	+4°C
Stop Solution	12 mL	+4°C
Sample Diluent NS	50 mL	+4°C
Anti-tag coated microplate (12 x 8 well strips)	96 Wells	+4°C
Plate Seal	1	+4°C

7. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Microplate reader capable of measuring absorbance at 450 or 600 nm.
- Method for determining protein concentration (BCA assay recommended).
- Deionized water.
- Multi- and single-channel pipettes.
- Tubes for standard dilution.
- Plate shaker for all incubation steps.
- Optional: Phenylmethylsulfonyl Fluoride (PMSF) (or other protease inhibitors).

8. Technical Hints

- Samples generating values higher than the highest standard should be further diluted in the appropriate sample dilution buffers.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions.
- Ensure plates are properly sealed or covered during incubation steps.
- Complete removal of all solutions and buffers during wash steps is necessary to minimize background.
- As a guide, typical ranges of sample concentration for commonly used sample types are shown below in Sample Preparation (section 11).
- All samples should be mixed thoroughly and gently.
- Avoid multiple freeze/thaw of samples.
- Incubate ELISA plates on a plate shaker during all incubation steps.
- When generating positive control samples, it is advisable to change pipette tips after each step.

- The provided Cell Extraction Enhancer Solution 50X may precipitate when stored at + 4°C. To dissolve, warm briefly at + 37°C and mix gently. The Cell Extraction Enhancer Solution 50X can be stored at room temperature to avoid precipitation.
- To avoid high background always add samples or standards to the well before the addition of the antibody cocktail.
- This kit is sold based on number of tests. A 'test' simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.

9. Reagent Preparation

- Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for 96 wells. The sample volumes below are sufficient for 48 wells (6 x 8-well strips); adjust volumes as needed for the number of strips in your experiment.
- Prepare only as much reagent as is needed on the day of the experiment. Capture and Detector Antibodies have only been tested for stability in the provided 10X formulations.

9.1 1X Cell Extraction Buffer PTR (For tissue extracts only):

Prepare 1X Cell Extraction Buffer PTR by diluting Cell Extraction Buffer PTR 5X and 50X Cell Extraction Enhancer Solution to 1X with deionized water. To make 10 mL 1X Cell Extraction Buffer PTR combine 7.8 mL deionized water, 2 mL Cell Extraction Buffer PTR 5X and 200 µL Cell Extraction Enhancer Solution 50X. Mix thoroughly and gently. If required protease inhibitors can be added.

Alternative – Enhancer may be added to 1X Cell Extraction Buffer PTR after extraction of cells or tissue. Refer to note in the Troubleshooting section.

9.2 1X Wash Buffer PT:

Prepare 1X Wash Buffer PT by diluting Wash Buffer PT 10X with deionized water. To make 50 mL 1X Wash Buffer PT combine 5 mL Wash Buffer PT 10X with 45 mL deionized water. Mix thoroughly and gently.

9.3 Antibody Cocktail:

Prepare Antibody Cocktail by diluting the capture and detector antibodies in Antibody Diluent 4BR. To make 3 mL of the Antibody Cocktail combine 300 µL 10X Capture Antibody and 300 µL 10X Detector Antibody with 2.4 mL Antibody Diluent 4BR. Mix thoroughly and gently.

10. Standard Preparation

- Always prepare a fresh set of standards for every use.
- Discard working standard dilutions after use as they do not store well.
- The following section describes the preparation of a standard curve for duplicate measurements (recommended).

10.1 **IMPORTANT:** If the protein standard vial has a volume identified on the label, reconstitute by adding that volume of dilutant indicated on the label. Alternatively, if the vial has a mass identified, reconstitute by adding 500 μL of dilutant.

10.2 For **serum, plasma, and tissue supernatant samples measurements**, reconstitute the CD93 protein standard in Sample Diluent NS.

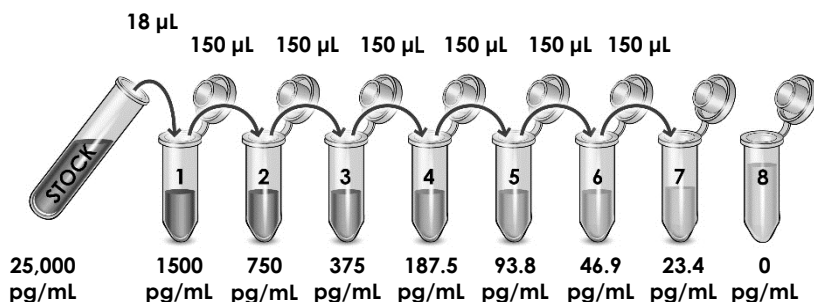
For **tissue homogenate extract samples measurements**, reconstitute the CD93 protein standard in 1X Cell Extraction Buffer PTR.

Hold at room temperature for 10 minutes and mix thoroughly and gently. This is the 25 ng/mL **Stock Standard Solution**.

10.3 Label eight tubes, Standards 1– 8.

10.4 Add 282 μL of appropriate diluent (see step 10.1) into tube number 1 and 150 μL of appropriate diluent into numbers 2-8.

10.5 Use the Stock Standard to prepare the following dilution series:



Add 180 μL of Stock Standard into Tube #1, then perform 1:2 serial dilutions for Tubes 2 - 7 as detailed above. Standard #8 contains no protein and is the Blank control.

11. Sample Preparation

Typical Sample Dynamic Range	
Sample Type	Range
Serum	1: 16,000 – 1: 1,000
Plasma – Citrate	1: 16,000 – 1: 500
Plasma – Heparin	1: 16,000 – 1: 1,000
Plasma – EDTA	1: 16,000 – 1: 500
Mouse lung supernatant	1: 6,400 – 1: 200
Mouse heart supernatant	1: 1600 – 1: 50
Mouse lung homogenate extract	0.1 – 3 µg/ml
Mouse heart homogenate extract	0.8 – 25 µg/ml
Mouse spleen homogenate extract	1.6 – 50 µg/ml

11.1 Serum:

Samples should be collected into a serum separator tube. After clot formation, centrifuge samples at 2,000 x g for 10 minutes and collect serum. Dilute samples into Sample Diluent NS and assay. Store un-diluted serum at -20°C or below. Avoid repeated freeze-thaw cycles.

11.2 Plasma:

Collect plasma using citrate, EDTA or heparin. Centrifuge samples at 2,000 x g for 10 minutes. Dilute samples into Sample Diluent NS and assay. Store un-diluted plasma samples at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

11.3 Tissue Culture Supernatants:

Centrifuge tissue culture media at 2,000 x g for 10 minutes to remove debris. Collect supernatants and assay. Or dilute samples into Sample Diluent NS and assay. Store un-diluted samples at -20°C or below. Avoid repeated freeze-thaw cycles.

11.4 Preparation of extracts from tissue homogenates:

- 11.4.1 Tissue lysates are typically prepared by homogenization of tissue that is first minced and thoroughly rinsed in PBS to remove blood (Dounce homogenizer recommended).
- 11.4.2 Homogenize 100 to 200 mg of wet tissue in 500 μ L – 1 mL of chilled 1X Cell Extraction Buffer PTR. For lower amounts of tissue adjust volumes accordingly.
- 11.4.3 Incubate on ice for 20 minutes.
- 11.4.4 Centrifuge at 18,000 x g for 20 minutes at 4°C.
- 11.4.5 Transfer the supernatants into clean tubes and discard the pellets.
- 11.4.6 Assay samples immediately or aliquot and store at -80°C. The sample protein concentration in the extract may be quantified using a protein assay.
- 11.4.7 Dilute samples to desired concentration in 1X Cell Extraction Buffer PTR.

12. Plate Preparation

- The 96 well plate strips included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents.
- Unused plate strips should be immediately returned to the foil pouch containing the desiccant pack, resealed and stored at 4°C.
- For each assay performed, a minimum of two wells must be used as the zero control.
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).
- Differences in well absorbance or "edge effects" have not been observed with this assay.

13. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature prior to use.
 - We recommend that you assay all standards, controls and samples in duplicate.
- 13.1 Prepare all reagents, working standards, and samples as directed in the previous sections.
 - 13.2 Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, reseal and return to 4°C storage.
 - 13.3 Add 50 μ L of all sample or standard to appropriate wells.
 - 13.4 Add 50 μ L of the Antibody Cocktail to each well.
 - 13.5 Seal the plate and incubate for 1 hour at room temperature on a plate shaker set to 400 rpm.
 - 13.6 Wash each well with 3 x 350 μ L 1X Wash Buffer PT. Wash by aspirating or decanting from wells then dispensing 350 μ L 1X Wash Buffer PT into each well. Wash Buffer PT should remain in wells for at least 10 seconds. Complete removal of liquid at each step is essential for good performance. After the last wash invert the plate and tap gently against clean paper towels to remove excess liquid.
 - 13.7 Add 100 μ L of TMB Development Solution to each well and incubate for 10 minutes in the dark on a plate shaker set to 400 rpm.

Given variability in laboratory environmental conditions, optimal incubation time may vary between 5 and 20 minutes.

Note: The addition of Stop Solution will change the color from blue to yellow and enhance the signal intensity about 3X. To avoid signal saturation, proceed to the next step before the high concentration of the standard reaches a blue color of O.D.600 equal to 1.0.
 - 13.8 Add 100 μ L of Stop Solution to each well. Shake plate on a plate shaker for 1 minute to mix. Record the OD at 450 nm. This is an endpoint reading.
 - 13.9 Alternative to 13.7 – 13.8: Instead of the endpoint reading at 450 nm, record the development of TMB Substrate kinetically. Immediately after addition of TMB Development Solution begin recording the blue color development with elapsed

time in the microplate reader prepared with the following settings:

Mode	Kinetic
Wavelength:	600 nm
Time:	up to 20 min
Interval:	20 sec - 1 min
Shaking:	Shake between readings

Δ **Note:** that an endpoint reading can also be recorded at the completion of the kinetic read by adding 100 µL Stop Solution to each well and recording the OD at 450 nm.

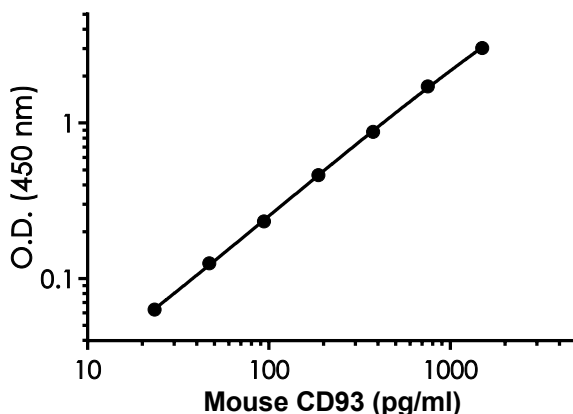
13.10 Analyze the data as described below.

14. Calculations

- 14.1 Calculate the average absorbance value for the blank control (zero) standards. Subtract the average blank control standard absorbance value from all other absorbance values.
 - 14.2 **Create a standard curve** by plotting the average blank control subtracted absorbance value for each standard concentration (y-axis) against the target protein concentration (x-axis) of the standard. Use graphing software to draw the best smooth curve through these points to construct the standard curve.
- Δ **Note:** Most microplate reader software or graphing software will plot these values and fit a curve to the data. A four-parameter curve fit (4PL) is often the best choice; however, other algorithms (e.g. linear, semi-log, log/log, 4-parameter logistic) can also be tested to determine if it provides a better curve fit to the standard values.
- 14.3 Determine the concentration of the target protein in the sample by interpolating the blank control subtracted **absorbance values against the standard curve**. Multiply the resulting value by the appropriate sample dilution factor, if used, to obtain the concentration of target protein in the sample.
 - 14.4 Samples generating absorbance values greater than that of the highest standard should be further diluted and reanalyzed. Similarly, samples which measure at an absorbance values less than that of the lowest standard should be retested in a less dilute form.

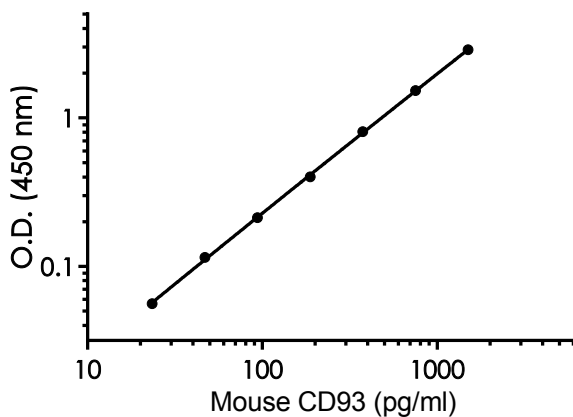
15. Typical Data

Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.



Standard Curve Measurements			
Concentration (pg/mL)	O.D 450 nm		Mean O.D
	1	2	
0	0.047	0.047	0.047
23.4	0.111	0.109	0.110
46.9	0.173	0.172	0.172
93.8	0.279	0.281	0.280
187.5	0.503	0.516	0.509
375	0.932	0.919	0.925
750	1.741	1.794	1.768
1500	3.047	3.121	3.084

Figure 1. Example of mouse CD93 standard curve in Sample Diluent NS. The CD93 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



Standard Curve Measurements			
Concentration (pg/mL)	O.D 450 nm		Mean O.D
	1	2	
0	0.047	0.047	0.047
23.4	0.101	0.105	0.103
46.9	0.160	0.164	0.162
93.8	0.257	0.262	0.259
187.5	0.442	0.453	0.447
375	0.870	0.837	0.853
750	1.556	1.595	1.576
1500	2.870	2.971	2.920

Figure 2. Example of mouse CD93 standard curve in 1X Cell Extraction Buffer PTR. The CD93 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

16. Calibration

This immunoassay is calibrated against a highly purified mouse CD93.

17. Typical Sample Values

SENSITIVITY –

The calculated minimal detectable dose (MDD) is 10.4 pg/mL in Sample Diluent NS and 3.4 pg/ml in 1X Cell Extraction Buffer PTR. The MDD was determined by calculating the mean of zero standard replicates (n=25) and adding 2 standard deviations then extrapolating the corresponding concentration.

RECOVERY –

Three concentrations of CD93 standard protein were spiked in duplicate to the indicated biological matrix to evaluate signal recovery in the working range of the assay.

Sample Type	Average % Recovery	Range (%)
Mouse Serum (1: 4,000)	102	98-107
Mouse Plasma – Citrate (1: 4,000)	103	99-110
Mouse Plasma – Heparin (1: 3,000)	108	101-113
Mouse Plasma – EDTA (1: 2,000)	101	98-106
Mouse lung supernatant (1: 1,000)	99	97-100
Mouse heart supernatant (1: 128)	103	97-112
Mouse lung homogenate extract (1 µg/ml)	108	102-117
Mouse heart homogenate extract (8 µg/ml)	94	86-101
Mouse spleen homogenate extract (10 µg/ml)	109	99-126%

Linearity of Dilution

Linearity of dilution is determined based on interpolated values from the standard curve. Linearity of dilution defines a sample concentration interval in which interpolated target concentrations are directly proportional to sample dilution.

Native CD93 was measured in the following biological samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

Dilution Factor	Interpolated value	1: 1,000 Mouse Serum	1: 500 Mouse Plasma (Citrate)	1: 1,000 Mouse Plasma (Heparin)	1: 500 Mouse Plasma (EDTA)
Undiluted	pg/mL	908	1264	795	940
	% Expected value	100	100	100	100
2	pg/mL	439	641	394	476
	% Expected value	97	101	99	101
4	pg/mL	213	318	206	247
	% Expected value	94	101	104	105
8	pg/mL	104	157	100	117
	% Expected value	92	99	101	100
16	pg/mL	53	79	49	56
	% Expected value	93	100	99	95

Native CD93 was measured in the following biological samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

Dilution Factor	Interpolated value	1: 200 Mouse Lung Supernatant	1: 50 Mouse Heart Supernatant
Undiluted	pg/mL	530	1019
	% Expected value	100	100
2	pg/mL	271	523
	% Expected value	102	103
4	pg/mL	126	257
	% Expected value	95	101
8	pg/mL	63	121
	% Expected value	95	95
16	pg/mL	31	62
	% Expected value	94	97

Native CD93 was measured in the following biological samples in a 2-fold dilution series. Sample dilutions are made in 1X Cell Extraction Buffer PTR.

Dilution Factor	Interpolated value	3 µg/ml Mouse Lung Homogenate Extract	25 µg/ml Mouse Heart Homogenate Extract	50 µg/ml Mouse Spleen Homogenate Extract
Undiluted	pg/mL	904	1133	1155
	% Expected value	100	100	100
2	pg/mL	435	544	582
	% Expected value	96	96	101
4	pg/mL	221	282	285
	% Expected value	98	100	99
8	pg/mL	110	137	135
	% Expected value	97	97	94
16	pg/mL	61	68	66
	% Expected value	108	96	91

PRECISION –

Mean coefficient of variations of interpolated values of CD93 from three concentrations of mouse serum within the working range of the assay.

	Intra- Assay	Inter- Assay
n =	5	3
CV (%)	4.2	5.7

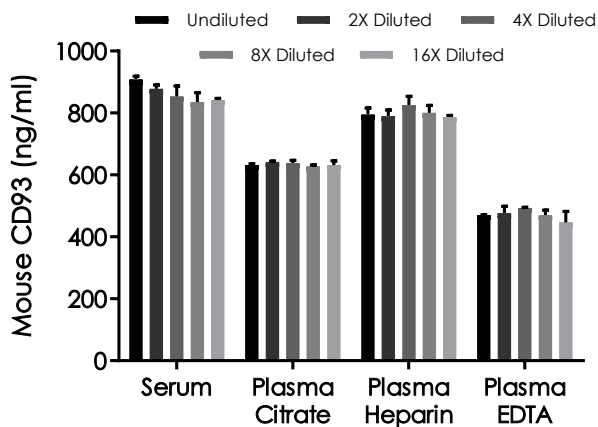


Figure 3. Interpolated concentrations of native CD93 in mouse serum and plasma samples. The concentrations of CD93 were measured in duplicates, interpolated from the CD93 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 1: 1,000, plasma (citrate) 1: 500, plasma (heparin) 1: 1,000, and plasma (EDTA) 1: 500. The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2). The mean CD93 concentration was determined to be 863 ng/mL in serum, 633 ng/mL in plasma (citrate), 799 ng/ml in plasma (heparin), and 468 ng/ml in plasma (EDTA).

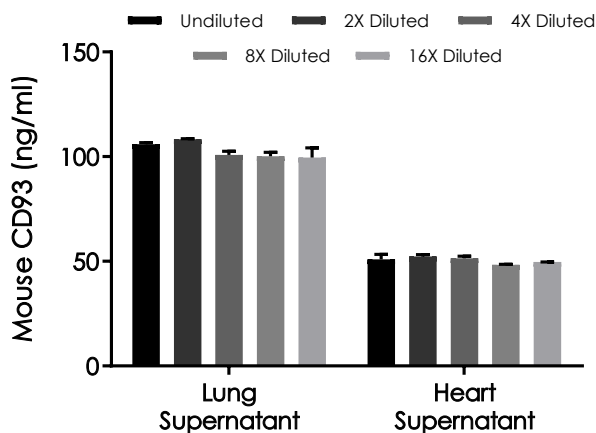


Figure 4. Interpolated concentrations of native CD93 in mouse tissue culture supernatant samples. The concentrations of CD93 were measured in duplicates, interpolated from the CD93 standard curves and corrected for sample dilution. Undiluted samples are as follows: mouse lung supernatant 1: 200 and mouse heart supernatant 2%. The interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean CD93 concentration was determined to be 103 ng/mL in mouse lung supernatant and 50 ng/ml in mouse heart supernatant.

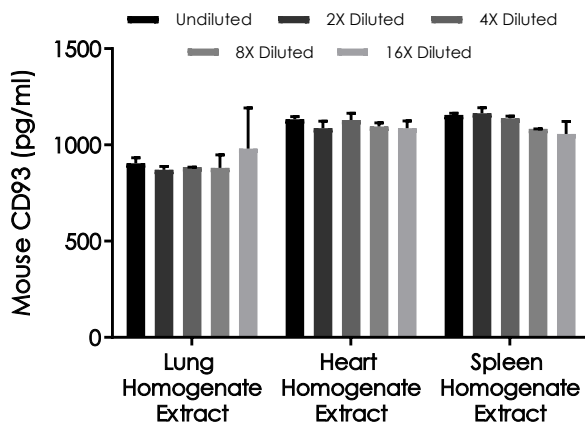


Figure 5. Interpolated concentrations of native CD93 in mouse lung, heart, and spleen homogenate extract samples based on a 3 µg/mL, 25 µg/mL, and 50 µg/mL extract load respectively. The concentrations of CD93 were measured in duplicate and interpolated from the CD93 standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CD93 concentration was determined to be 898 pg/mL in lung homogenate extract, 1116 pg/ml in heart homogenate extract, and 1133 ng/mL in spleen homogenate extract samples.

18. Assay Specificity

This kit recognizes both native and recombinant mouse CD93 protein in serum, plasma, and cell culture supernatant, cell and tissue extract samples only.

Cell culture supernatant and cell culture extract samples have not been tested with this kit.

CROSS REACTIVITY

Recombinant human CD93 was prepared at 50 ng/mL and 1.5 ng/mL and assayed for cross reactivity. No cross-reactivity was observed.

INTERFERENCE

Recombinant mouse MBL-2 was prepared at 50 ng/mL and 1.5 ng/mL and tested for interference. No interference with was observed.

19. Species Reactivity

This kit recognizes mouse CD93 protein.

Other species reactivity was determined by measuring 1: 2,000 human, rat, and bovine serum samples, interpolating the mouse CD93 protein concentrations from the mouse standard curve, and expressing the interpolated concentrations as a percentage of the mouse CD93 protein concentration in mouse serum assayed at the same dilution.

No cross-reactivity was observed.

Please contact our Technical Support team for more information.

20. Troubleshooting

Problem	Reason	Solution
Difficulty pipetting lysate; viscous lysate.	Genomic DNA solubilized	Prepare 1X Cell Extraction Buffer PTR (without enhancer). Add enhancer to lysate after extraction.
Poor standard curve	Inaccurate Pipetting	Check pipettes
	Improper standard dilution	Prior to opening, briefly spin the stock standard tube and dissolve the powder thoroughly by gentle mixing
Low Signal	Incubation times too brief	Ensure sufficient incubation times; increase to 2 or 3 hour standard/sample incubation
	Inadequate reagent volumes or improper dilution	Check pipettes and ensure correct preparation
	Incubation times with TMB too brief	Ensure sufficient incubation time until blue color develops prior addition of Stop solution
Large CV	Plate is insufficiently washed	Review manual for proper wash technique. If using a plate washer, check all ports for obstructions.
	Contaminated wash buffer	Prepare fresh wash buffer
Low sensitivity	Improper storage of the ELISA kit	Store your reconstituted standards at -80°C, all other assay components 4°C. Keep TMB Development Solution protected from light.
Precipitate in Diluent	Precipitation and/or coagulation of components within the Diluent.	Precipitate can be removed by gently warming the Diluent to 37°C.

Technical Support

Copyright © 2023 Abcam, All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)