



ab203365 - Osteoprotegerin (TNFRSF11B) Mouse SimpleStep ELISA[®] Kit

Instructions for Use

For the quantitative measurement of Osteoprotegerin (TNFRSF11B) in mouse serum, plasma and cell culture supernatants samples.

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

Table of Contents

INTRODUCTION

1. BACKGROUND	2
2. ASSAY SUMMARY	3

GENERAL INFORMATION

3. PRECAUTIONS	4
4. STORAGE AND STABILITY	4
5. MATERIALS SUPPLIED	4
6. MATERIALS REQUIRED, NOT SUPPLIED	4
7. LIMITATIONS	5
8. TECHNICAL HINTS	5

ASSAY PREPARATION

9. REAGENT PREPARATION	7
10. STANDARD PREPARATION	8
11. SAMPLE PREPARATION	10
12. PLATE PREPARATION	11

ASSAY PROCEDURE

13. ASSAY PROCEDURE	12
---------------------	----

DATA ANALYSIS

14. CALCULATIONS	14
15. TYPICAL DATA	15
16. TYPICAL SAMPLE VALUES	17
17. ASSAY SPECIFICITY	21
18. SPECIES REACTIVITY	22

RESOURCES

19. TROUBLESHOOTING	23
20. NOTES	24

1. BACKGROUND

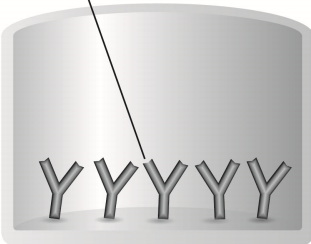
Abcam's Osteoprotegerin (TNFRSF11B) *in vitro* SimpleStep ELISA® (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Osteoprotegerin protein in mouse serum, plasma and cell culture supernatants samples.

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.

Osteoprotegerin acts as decoy receptor for TNFSF11/RANKL and thereby neutralizes its function in osteoclastogenesis. Osteoprotegerin inhibits the activation of osteoclasts and promotes osteoclast apoptosis *in vitro*. Bone homeostasis seems to depend on the local ratio between TNFSF11 and TNFRSF11B. Osteoprotegerin may also play a role in preventing arterial calcification. Osteoprotegerin may act as decoy receptor for TNFSF10/TRAIL and protect against apoptosis. TNFSF10/TRAIL binding blocks the inhibition of osteoclastogenesis.

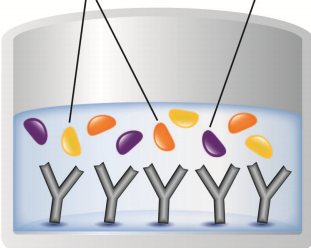
2. ASSAY SUMMARY

Immobilization Antibody



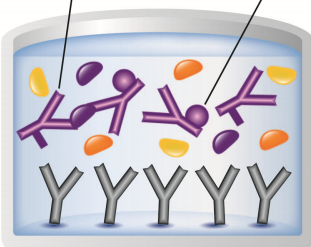
Remove appropriate number of antibody coated well strips. Equilibrate all reagents to room temperature. Prepare all reagents, samples, and standards as instructed.

Matrix Proteins Target Analyte



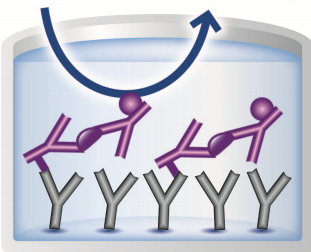
Add standard or sample to appropriate wells.

Capture Antibody Detector Antibody



Add Antibody Cocktail to all wells. Incubate at room temperature.

Substrate Color Development



Aspirate and wash each well. Add TMB Development Solution to each well and incubate. Add Stop Solution at a defined endpoint. Alternatively, record color development kinetically after TMB substrate addition.

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. Modifications to the kit components or procedures may result in loss of performance.

4. STORAGE AND STABILITY

Store kit at 2-8°C immediately upon receipt.

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Reagent and Standard Preparation sections.

5. MATERIALS SUPPLIED

Item	Amount	Storage Condition (Before Preparation)
10X Mouse Osteoprotegerin Capture Antibody	600 µL	+2-8°C
10X Mouse Osteoprotegerin Detector Antibody	600 µL	+2-8°C
Mouse Osteoprotegerin Lyophilized Recombinant Protein	2 Vials	+2-8°C
Antibody Diluent 4BI	6 mL	+2-8°C
10X Wash Buffer PT	20 mL	+2-8°C
TMB Development Solution	12 mL	+2-8°C
Stop Solution	12 mL	+2-8°C
Sample Diluent NS	50 mL	+2-8°C
Sample Diluent 50BS	20 mL	+2-8°C
Pre-Coated 96 Well Microplate (12 x 8 well strips)	96 Wells	+2-8°C
Plate Seal	1	+2-8°C

6. MATERIALS REQUIRED, NOT SUPPLIED

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

- Microplate reader capable of measuring absorbance at 450 or 600 nm.
- Method for determining protein concentration (BCA assay recommended).
- Deionized water.
- PBS (1.4 mM KH₂PO₄, 8 mM Na₂HPO₄, 140 mM NaCl, 2.7 mM KCl, pH 7.4).
- Multi- and single-channel pipettes.
- Tubes for standard dilution.
- Plate shaker for all incubation steps.
- Phenylmethylsulfonyl Fluoride (PMSF) (or other protease inhibitors).

7. 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.

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.
- **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 Wash Buffer PT**

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

9.2 **Antibody Cocktail**

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

10. STANDARD PREPARATION

Prepare serially diluted standards immediately prior to use. Always prepare a fresh set of positive controls for every use.

The following section describes the preparation of a standard curve for duplicate measurements (recommended).

IMPORTANT: If the protein standard vial has a volume identified on the label, reconstitute the Osteoprotegerin standard by adding that volume of Diluent indicated on the label. Alternatively, if the vial has a mass identified, reconstitute the Osteoprotegerin standard by adding 200 μ L Diluent. Hold at room temperature for 10 minutes and mix gently. This is the 16,000 pg/mL **Stock Standard** Solution.

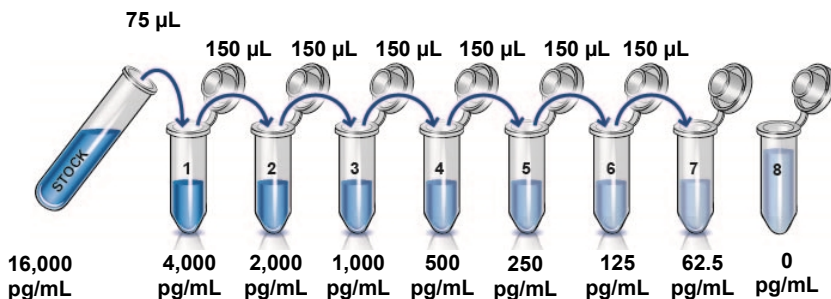
10.1 For **cell culture supernatant samples** follow this section to prepare the standard.

10.1.1 Reconstitute the Osteoprotegerin standard by adding water by pipette.

10.1.2 Label eight tubes, Standards 1–8.

10.1.3 Add 225 μ L Sample Diluent NS into tube number 1 and 150 μ L of Sample Diluent NS into numbers 2–8.

10.1.4 Use the Stock Standard to prepare the following dilution series. Standard #8 contains no protein and is the Blank control:



ASSAY PROCEDURE

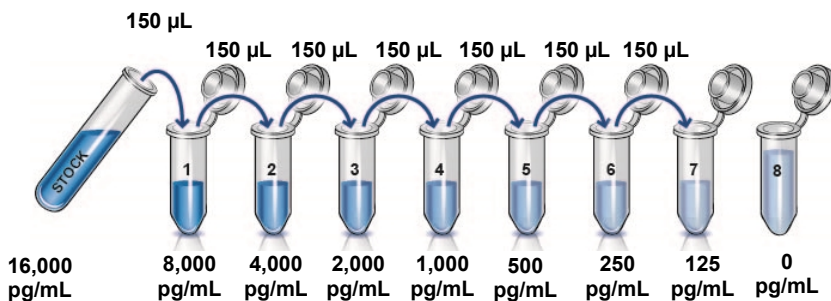
10.2 For **serum and plasma (heparin) samples** follow this section.

10.2.1 Reconstitute the Osteoprotegerin standard by adding Sample Diluent 50BS by pipette.

10.2.2 Label eight tubes, Standards 1– 8.

10.2.3 Add 150 μ L Sample Diluent 50BS into tube number 1 and 150 μ L of Sample Diluent 50BS into numbers 2-8.

10.2.4 Use the Stock Standard to prepare the following dilution series. Standard #8 contains no protein and is the Blank control:



11. SAMPLE PREPARATION

TYPICAL SAMPLE DYNAMIC RANGE	
Sample Type	Range
Mouse lung day 5 cell culture supernatant	Neat - 25%
Mouse serum	Neat - 25%
Mouse plasma (heparin)	Neat - 12.5%

11.1 Plasma

Collect plasma using heparin. Centrifuge samples at 2,000 x g for 10 minutes. Plasma (heparin) samples can be assayed without dilution. If needed, dilute samples into Sample Diluent 50BS and assay. Store un-diluted plasma samples at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

11.2 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. Serum samples can be assayed without dilution. If needed, dilute samples into Sample Diluent 50BS and assay. Store un-diluted serum at -20°C or below. Avoid repeated freeze-thaw cycles.

11.3 Cell Culture Supernatants

Centrifuge cell culture media at 2,000 x g for 10 minutes to remove debris. Collect supernatants. Cell culture supernatant samples can be assayed without dilution. If needed, dilute samples into Sample Diluent NS and assay. Store samples at -20°C or below. Avoid repeated freeze-thaw cycles.

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.**
- **It is recommended to 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 µL of all sample or standard to appropriate wells.
- 13.4 Add 50 µL 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 µL 1X Wash Buffer PT. Wash by aspirating or decanting from wells then dispensing 350 µL 1X Wash Buffer PT into each well. Complete removal of liquid at each step is essential for good performance. After the last wash invert the plate and blot it against clean paper towels to remove excess liquid.
- 13.7 Add 100 µL 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 µL 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.

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

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

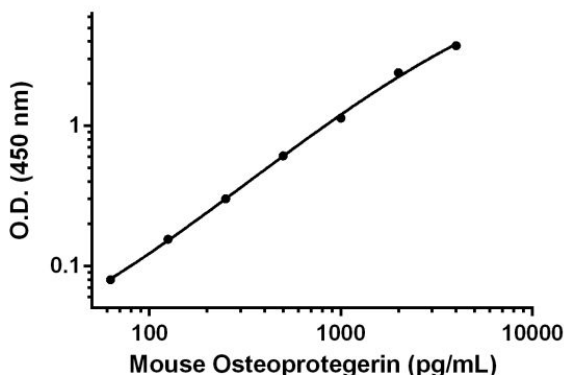
13.9 Analyze the data as described below.

14. CALCULATIONS

Subtract the average zero standard absorbance value from all measurements to determine the background-corrected absorbance value. Determine the average value of the duplicate background-subtracted measurements of the positive control dilutions and plot these values against their concentrations. Draw the best smooth curve through these points to construct a standard curve. Software supplied with microplate readers have or graphing software can plot these values and curve fit. A four parameter algorithm (4PL) usually provides the best fit, though other equations can be examined to determine the most accurate (e.g. linear, semi-log, log/log, 4 parameter logistic). The concentrations of unknown samples can be interpolated from the curve using the background-corrected absorbance values and the dilution factor, if used. Absorbance values of samples which are greater than that of the highest standard should be diluted until the absorbance value falls within the linear part of the curve. The concentration of the sample can then be determined by applying the dilution factor.

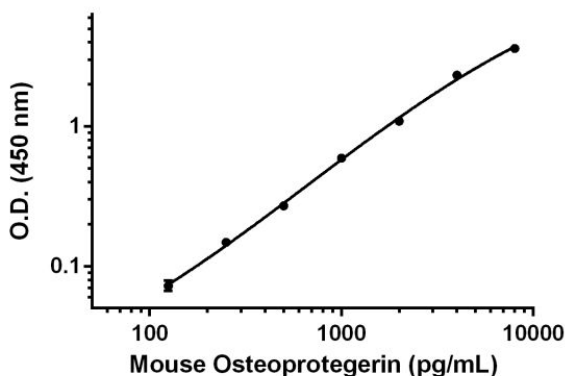
15. TYPICAL DATA

TYPICAL STANDARD CURVE – Data provided for **demonstration purposes only**. A new standard curve must be generated for each assay performed.



Standard Curve Measurements			
Conc. (pg/mL)	O.D. 450 nm		Mean O.D.
	1	2	
0	0.051	0.053	0.052
62.5	0.131	0.133	0.132
125	0.203	0.211	0.207
250	0.348	0.359	0.354
500	0.637	0.688	0.663
1,000	1.156	1.222	1.190
2,000	2.467	2.431	2.449
4,000	3.806	3.772	3.789

Figure 1. Example of Osteoprotegerin standard curve for prepared in Sample Diluent NS measurements. The Osteoprotegerin 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			
Conc. (pg/mL)	O.D. 450 nm		Mean O.D.
	1	2	
0	0.050	0.047	0.049
125	0.126	0.117	0.122
250	0.197	0.196	0.197
500	0.319	0.320	0.320
1,000	0.665	0.619	0.642
2,000	1.188	1.095	1.142
4,000	2.478	2.273	2.376
8,000	3.681	3.650	3.666

Figure 2. Example of Osteoprotegerin standard curve for prepared in Sample Diluent 50BS measurements. The Osteoprotegerin 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. TYPICAL SAMPLE VALUES

SENSITIVITY –

The calculated minimal detectable dose (MDD) is determined by calculating the mean of zero standard replicates and adding 2 standard deviations then extrapolating the corresponding concentrations. The MDD is dependent on the Sample Diluent buffer used:

Sample Diluent Buffer	n=	Minimal Detectable Dose
Sample Diluent NS	22	13 pg/mL
Sample Diluent 50BS	25	5 pg/mL

RECOVERY –

Three concentrations of Mouse Recombinant Osteoprotegerin 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 (%)
100% Cell Culture Media	104.9	102.1 – 106.3
100% Mouse serum	96.1	90.0 – 101.3
100% Mouse plasma (heparin)	123.4	117.9 – 132.8

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 Osteoprotegerin was measured in the following biological samples in a 2-fold dilution series. Cell culture supernatant sample dilutions were made in Sample Diluent NS. Serum and plasma (heparin) sample dilutions were made in Sample Diluent 50BS.

DATA ANALYSIS

Dilution Factor	Interpolated value	100% Mouse Lung Day 5 Supern.	100% Mouse Serum	100% Mouse Plasma (Heparin)
Undiluted	pg/mL	374.98	1,136.03	1,062.78
	% Expected value	100	100	100
2	pg/mL	206.59	550.65	600.72
	% Expected value	110.2	96.9	113.1
4	pg/mL	91.08	243.97	300.10
	% Expected value	97.2	85.9	113.0
8	pg/mL	NA	NA	125.04
	% Expected value	NA	NA	94.1

NA – Not Analyzed, measured O.D. values were lower than the O.D. value of the lowest standard

Mouse recombinant Osteoprotegerin was spiked into the following biological samples and diluted in a 2-fold dilution series. Cell culture media (10F RPMI1640) samples were diluted in Sample Diluent NS. Mouse serum and plasma (heparin) samples were diluted in Sample Diluent 50BS.

Dilution Factor	Interpolated value	100% Cell Culture Media	100% Mouse Serum	100% Mouse Plasma (Heparin)
Undiluted	pg/mL	2,550.07	6,991.94	6,916.99
	% Expected value	100	100	100
2	pg/mL	1,264.21	3,261.05	3,636.62
	% Expected value	99.2	93.3	105.2
4	pg/mL	656.13	1,422.70	1,445.43
	% Expected value	102.9	81.4	83.6
8	pg/mL	335.77	776.41	757.19
	% Expected value	105.3	88.8	87.6
16	pg/mL	155.19	395.50	368.26
	% Expected value	97.4	90.5	85.2
32	pg/mL	68.85	206.41	173.95
	% Expected value	86.4	94.5	80.5

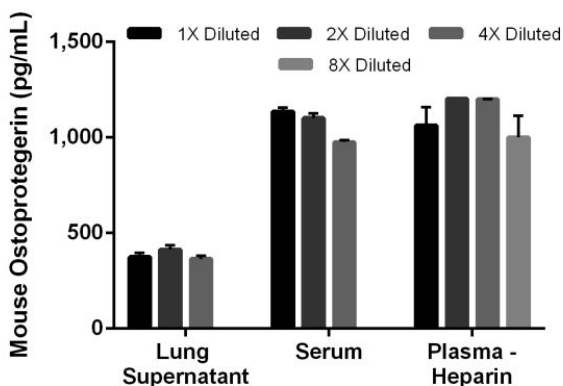


Figure 3. Interpolated concentrations of Osteoprotegerin in mouse lung day 5 supernatant, serum and plasma (heparin) samples. The concentrations of Osteoprotegerin were measured in duplicates, interpolated from the Osteoprotegerin standard curve and corrected for sample dilution. Note that 1X Diluted samples were neat. The interpolated, dilution factor-corrected values are plotted in pg of Osteoprotegerin per mL of neat sample (mean $\pm$ SD, n=2).

Serum and Plasma. Pooled female mouse serum and plasma samples were measured in three different dilutions in duplicates for Osteoprotegerin concentrations using this kit. Means of interpolated sample values adjusted to sample dilution are shown in pg of Osteoprotegerin per mL of serum/plasma in the table below.

Sample Type	Mean (pg/mL)	Range (pg/mL)
Mouse Serum	1,071	976 – 1,136
Mouse Plasma - Heparin	1,155	1,063 – 1,201

PRECISION –

Mean coefficient of variations of interpolated values from 3 concentrations of mouse serum within the working range of the assay.

	Intra-Assay	Inter-Assay
n=	5	3
CV (%)	4.9	8.1

17. ASSAY SPECIFICITY

This kit recognizes both native and recombinant mouse Osteoprotegerin protein in serum, plasma (heparin) and cell culture supernatant samples only. Plasma (citrate and EDTA) samples are not compatible with this kit.

Urine, milk, saliva, and cell and tissue extract samples have not been tested with this kit.

CROSS REACTIVITY

Recombinant mouse TNF Receptor I and TNF Receptor II proteins were prepared at 8,000, 4,000, 2,000 and 1,000 pg/mL and assayed for cross reactivity. No cross-reactivity was observed.

18. SPECIES REACTIVITY

This kit recognizes mouse Osteoprotegerin protein.

Other species reactivity was determined by measuring neat serum samples of various species, interpolating the protein concentrations from the mouse standard curve, and expressing the interpolated concentrations as a percentage of the protein concentration in mouse serum assayed at the same dilution.

Reactivity < 3% was determined for the following species:

- Rat
- Rabbit
- Dog
- Goat
- Pig
- Cow
- Human

Please contact our Scientific Support team for more information

19. TROUBLESHOOTING

Problem	Cause	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.

20. NOTES

Technical Support

Copyright © 2021 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)