

Version 2 Last updated 18 November 2021

ab264619 Rat Beta-2- Microglobulin SimpleStep ELISA[®] Kit

For the quantitative measurement of Beta-2-Microglobulin in rat serum, plasma, urine, and cell culture supernatant 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	12
13. Assay Procedure	13
14. Calculations	15
15. Typical Data	16
16. Typical Sample Values	17
17. Assay Specificity	23
18. Species Reactivity	23
19. Troubleshooting	24
20. Notes	25
Technical Support	26

1. Overview

Beta-2-Microglobulin *in vitro* SimpleStep ELISA® (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Beta-2-Microglobulin protein in rat serum, plasma, urine, and cell culture supernatant 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.

Beta-2-Microglobulin is the light chain of the class I major histocompatibility complex (MHC); it forms heterodimers with classical (HLA-A, -B, -C) and non-classical (HLA-E, -F, -G) class I MHC heavy chains and with class I non-MHC heavy chains (CD1). Beta-2-Microglobulin is involved in the presentation of peptide antigens to the immune system. Beta-2-Microglobulin can dissociate from the MHC and circulate in complex with alpha 2-Macroglobulin.

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
Rat Beta-2-Microglobulin Capture Antibody 10X	600 µL	+4°C
Rat Beta-2-Microglobulin Detector Antibody 10X	600 µL	+4°C
Rat Beta-2-Microglobulin Lyophilized Recombinant Protein	2 Vials	+4°C
Antibody Diluent 5BR	6 mL	+4°C
Cell Extraction Enhancer Solution 50X	1 mL	+4°C
Denaturant	500 µL	+4°C
Sample Diluent NS	50 mL	+4°C
Wash Buffer PT 10X	20 mL	+4°C
TMB Development Solution	12 mL	+4°C
Stop Solution	12 mL	+4°C
SimpleStep Pre-Coated 96-Well Microplate	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 Sample Diluent NS + Enhancer:

Prepare Sample Diluent NS + Enhancer by diluting Cell Extraction Enhancer Solution 50X to 1X with Sample Diluent NS. To make 10 mL Sample Diluent NS + Enhancer combine 9.8 mL Sample Diluent NS and 200 μ L Cell Extraction Enhancer Solution 50X. Mix thoroughly and gently.

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 5BR. 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 5BR. Mix thoroughly and gently.

10. Standard Preparation

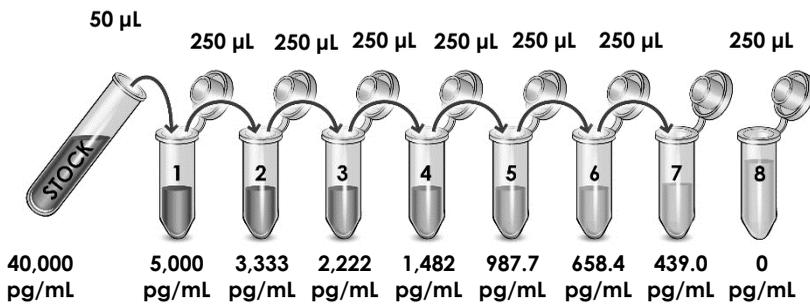
- Always prepare a fresh set of standards for every use.
- Discard working standard dilutions after use as they do not store well.
- The following section describes the preparation of a standard curve for duplicate measurements (recommended).

10.1 IMPORTANT: If the protein standard vial has a volume identified on the label, reconstitute the Beta-2-Microglobulin by adding that volume of Sample Diluent NS + Enhancer indicated on the label. Alternatively, if the vial has a mass identified, reconstitute the Beta-2-Microglobulin standard by adding 500 μ L of Sample Diluent NS + Enhancer. Hold at room temperature for 10 minutes and mix gently. This is the 40,000 pg/mL **Stock Standard** Solution.

10.1.1 Label eight tubes, Standards 1– 8.

10.1.2 Add 350 μ L Sample Diluent NS + Enhancer into tube number 1 and 125 μ L of Sample Diluent NS + Enhancer into numbers 2- 8.

10.1.3 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
Serum	1:22,781 - 1:3,000
Plasma - Citrate	1:11,391 - 1:1,500
Plasma - Heparin	1:17,086 - 1:2,250
Plasma - EDTA	1:11,391 - 1:2,250
Urine	1:759 - 1:150
Lung Cell Culture Supernatant	1:759 - 1:100

Note: The following dilution schemes have built-in required sample denaturation and 40-fold dilution of the Denaturant. As an example, the table below demonstrates the steps suggested to generate a final sample dilution. Final dilutions are suggestions based on observed sample reactivity and can be adjusted as needed.

For serum and plasma samples:

Tube #	Sample to Dilute	Volume to Dilute (μL)	Volume of Sample Diluent NS + Enhancer (μL)	Volume of Denaturant (μL)	Volume of Sample Diluent NS (μL)	Starting Dilution	Final Dilution
1	Neat Serum or Plasma	44.5	0	5.5	0	Neat	1:1.25
2	Tube #1	5	0	0	195	1:1.25	1:44.9

3	Tube #2	6	394	0	0	1:44.9	1:3,000
---	------------	---	-----	---	---	--------	---------

For urine samples:

Tube #	Sample to Dilute	Volume to Dilute (µL)	Volume of Sample Diluent NS + Enhancer (µL)	Volume of Denaturant (µL)	Volume of Sample Diluent NS (µL)	Starting Dilution	Final Dilution
1	Neat Urine	44.5	0	5.5	0	Neat	1:1.25
2	Tube #1	5	0	0	195	1:1.25	1:44.9
3	Tube #2	60	140	0	0	1:44.9	1:150

For cell culture supernatant samples:

Tube #	Sample to Dilute	Volume to Dilute (µL)	Volume of Sample Diluent NS + Enhancer (µL)	Volume of Denaturant (µL)	Volume of Sample Diluent NS (µL)	Starting Dilution	Final Dilution
1	Neat cell culture supernatant	44.5	0	5.5	0	Neat	1:1.25
2	Tube #1	5	0	0	195	1:1.25	1:44.9
3	Tube #2	90	110	0	0	1:44.9	1:100

11.1 Plasma:

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

- 11.1.1 To 44.5 volume parts of plasma add 5.5 volume part of Denaturant. Mix thoroughly and gently.
- 11.1.2 Incubate samples at room temperature for 10 minutes.
- 11.1.3 Dilute samples 40-fold in Sample Diluent NS. Mix thoroughly and gently.
- 11.1.4 Dilute samples to a final concentration of at least 1:2,250 in Sample Diluent NS + Enhancer and assay.

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. Store un-diluted serum at -20°C or below. Avoid repeated freeze-thaw cycles.

- 11.2.1 To 44.5 volume parts of serum add 5.5 volume part of Denaturant. Mix thoroughly and gently.
- 11.2.2 Incubate samples at room temperature for 10 minutes.
- 11.2.3 Dilute samples 40-fold in Sample Diluent NS. Mix thoroughly and gently.
- 11.2.4 Dilute samples to a final concentration of at least 1:3,000 in Sample Diluent NS + Enhancer and assay.

11.3 Cell Culture Supernatants:

Centrifuge cell culture media at 2,000 x g for 10 minutes to remove debris. Collect supernatants. Store un-diluted samples at -20°C or below. Avoid repeated freeze-thaw cycles.

- 11.3.1 To 44.5 volume parts of serum add 5.5 volume part of Denaturant. Mix thoroughly and gently.
- 11.3.2 Incubate samples at room temperature for 10 minutes.
- 11.3.3 Dilute samples 40-fold in Sample Diluent NS. Mix thoroughly and gently.
- 11.3.4 Assay or dilute samples in Sample Diluent NS + Enhancer and assay.

11.4 Urine:

Centrifuge urine at 2,000 x g for 10 minutes to remove debris. Store un-diluted urine samples at -20°C or below. Avoid repeated freeze-thaw cycles.

- 11.4.1 To 44.5 volume parts of serum add 5.5 volume part of Denaturant. Mix thoroughly and gently.
- 11.4.2 Incubate samples at room temperature for 10 minutes.
- 11.4.3 Dilute samples 40-fold in Sample Diluent NS. Mix thoroughly and gently.
- 11.4.4 Dilute samples to a final concentration of at least 1:150 in Sample Diluent NS + Enhancer and assay.

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.

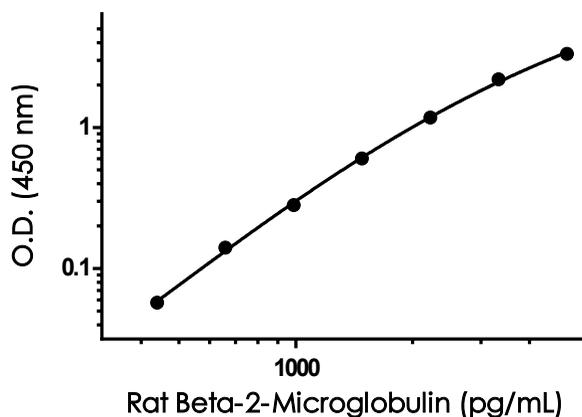
13.10 Analyze the data as described below.

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	0.070	0.074	0.072
439.0	0.127	0.131	0.129
658.4	0.212	0.214	0.213
987.7	0.358	0.349	0.354
1,482	0.680	0.672	0.676
2,222	1.254	1.246	1.250
3,333	2.237	2.312	2.274
5,000	3.408	3.402	3.405

Figure 1. Example of rat Beta-2-Microglobulin standard curve in Sample Diluent NS + Enhancer. The Beta-2-Microglobulin 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 268 pg/mL. The MDD was determined by calculating the mean of zero standard replicates (n=15) and adding 2 standard deviations then extrapolating the corresponding concentration.

RECOVERY –

Three concentrations of Beta-2-Microglobulin recombinant 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 (%)
1:10,000 Serum	115	111 - 118
1:10,000 Plasma - Citrate	109	107 - 111
1:10,000 Plasma - Heparin	111	106 - 114
1:10,000 Plasma - EDTA	118	117 - 119
1:250 Urine	119	116 - 123
1:333 Lung Cell Culture Supernatant	115	110 - 123
1:50 Cell Culture Media	99	93 - 104

*Media is RPMI 1640 containing 10% fetal bovine serum.

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 Beta-2-Microglobulin was measured in the following biological samples in a 1.5-fold dilution series. Sample dilutions are made in Sample Diluent NS + Enhancer.

Dilution Factor	Interpolated value	1:3,000 Rat Serum	1:1,500 Rat Plasma (Citrate)	1:2,250 Rat Plasma (Heparin)	1:2,250 Rat Plasma (EDTA)
Undiluted	pg/mL	4,451	4,745	4,421	3,472
	% Expected value	100	100	100	100
1.5	pg/mL	2,869	3,298	2,673	2,175
	% Expected value	97	104	91	94
2.25	pg/mL	1,785	2,136	1,698	1,431
	% Expected value	90	101	86	93
3.38	pg/mL	1,147	1,348	1,073	916.6
	% Expected value	87	96	82	89
5.06	pg/mL	738.1	833.2	688.4	563.5
	% Expected value	84	89	79	82

Native Beta-2-Microglobulin was measured in the following biological samples in a 1.5-fold dilution series. Sample dilutions are made in Sample Diluent NS + Enhancer.

Dilution Factor	Interpolated value	1:150 Rat Urine	1:100 Rat Lung Cell Culture Supernatant
Undiluted	pg/mL	3,418	4,395
	% Expected value	103	100
1.5	pg/mL	2,255	3,074
	% Expected value	102	105
2.25	pg/mL	1,437	1,857
	% Expected value	98	95
3.38	pg/mL	943.6	1,278
	% Expected value	96	98
5.06	pg/mL	625.6	832.4
	% Expected value	96	96

PRECISION –

Mean coefficient of variations of interpolated values of Beta-2-Microglobulin from two concentrations of rat serum within the working range of the assay.

	Intra- Assay	Inter- Assay
n =	8	3
CV (%)	3.2	2.1

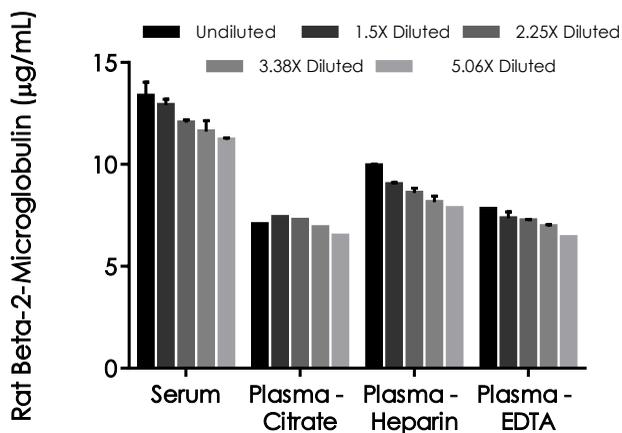


Figure 2. Interpolated concentrations of native Beta-2-Microglobulin in rat serum and plasma samples. The concentrations of Beta-2-Microglobulin were measured in duplicates, interpolated from the Beta-2-Microglobulin standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 1:3,000, plasma (citrate) 1:1,500, plasma (heparin) 1:2,250, and plasma (EDTA) 1:2,250. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Beta-2-Microglobulin concentration was determined to be 12.23 µg/mL in neat serum, 7.025 µg/mL in neat plasma (citrate), 8.710 µg/mL in neat plasma (heparin), and 7.155 µg/mL in neat plasma (EDTA).

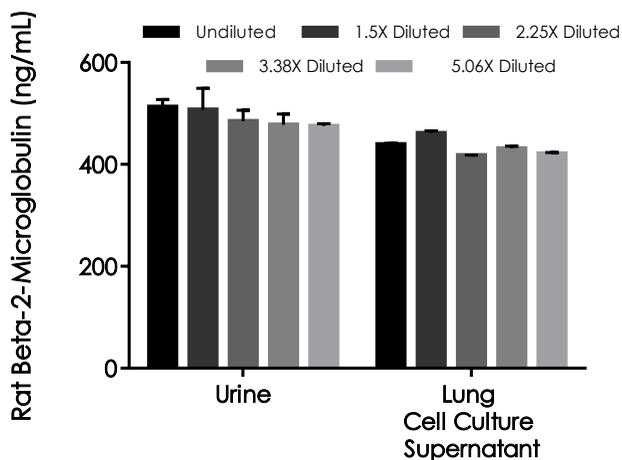


Figure 3. Interpolated concentrations of native Beta-2-Microglobulin in rat urine and cell culture supernatant samples. The concentrations of Beta-2-Microglobulin were measured in duplicates, interpolated from the Beta-2-Microglobulin standard curves and corrected for sample dilution. Undiluted samples are as follows: urine 1:150 and lung cell culture supernatant 1:100. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Beta-2-Microglobulin concentration was determined to be 491.5 ng/mL in neat rat urine and 434.3 ng/mL in neat rat lung cell culture supernatant.

17. Assay Specificity

This kit recognizes both native and recombinant rat Beta-2-Microglobulin protein in serum, plasma, and cell culture supernatant, and urine samples only.

Cell and tissue extract samples have not been tested with this kit.

18. Species Reactivity

This kit recognizes rat Beta-2-Microglobulin protein.

Other species reactivity was determined by measuring 1:5,000 serum samples of various species, interpolating the protein concentrations from the rat standard curve, and expressing the interpolated concentrations as a percentage of the protein concentration in rat serum assayed at the same dilution.

Reactivity < 3% was determined for the following species:

- Human
- Monkey
- Cow

This kit is not compatible with mouse samples.

Please contact our Technical Support team for more information.

19. 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.

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)