



ab199081 – IL-17A
Mouse SimpleStep
ELISA[®] Kit

Instructions for Use

For the quantitative measurement of IL-17A in mouse serum, plasma (citrate) and cell culture supernatant 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 5
- 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 21

RESOURCES

- 19. TROUBLESHOOTING 23
- 20. NOTES 24

1. BACKGROUND

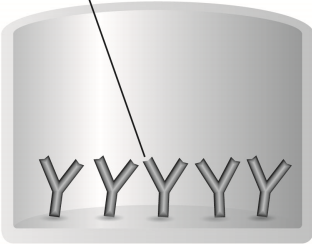
IL-17A (Interleukin-17a) mouse *in vitro* SimpleStep ELISA® (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-17A protein in mouse serum, plasma (citrate) and cell culture supernatant samples.

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.

IL-17A is a pro-inflammatory cytokine that is secreted by a subset of activated T cells. It is a disulfide-linked homodimer with both glycosylated and non-glycosylated forms. IL-17A induces stromal cells to produce pro-inflammatory and hematopoietic cytokines, and also enhances the surface expression of ICAM1/intracellular adhesion molecule 1 in fibroblasts.

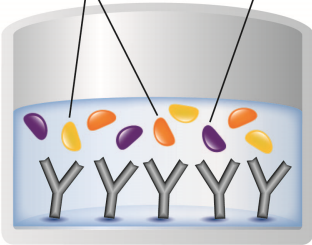
2. ASSAY SUMMARY

Immobilization Antibody



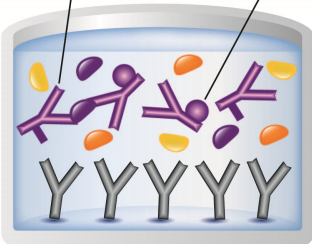
Remove appropriate number of antibody coated well strips. Equilibrate all reagents to room temperature. Prepare all reagents, samples, and standards as instructed.

Matrix Proteins Target Analyte



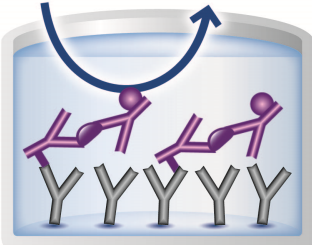
Add standard or sample to appropriate wells.

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 IL-17A Capture Antibody	600 µL	+2-8°C
10X Mouse IL-17A Detector Antibody	600 µL	+2-8°C
Mouse IL-17A Lyophilized Recombinant Protein	2 Vials	+2-8°C
Antibody Diluent CPI2	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 25BP	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

Note: Antibody Diluent CPI2- This buffer has been reformulated to enhance stability after freeze-thaw cycles while producing data equivalent to the original formulation of antibody diluent CPI previously used in this kit.

While we run stock down, you may receive kits containing antibody diluent CPI. This does not affect the way you should use the kit.

If you have any questions please contact Abcam Scientific Support.

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 KH_2PO_4 , 8 mM Na_2HPO_4 , 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
- Optional: 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 CPI2. 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 CPI2. 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 table 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 IL-17A standard by adding that volume of Diluent indicated on the label. Alternatively, if the vial has a mass identified, reconstitute the IL-17A standard by adding 200 μL Diluent. Hold at room temperature for 10 minutes and mix gently. This is the 1,600 pg/mL **Stock Standard** Solution.

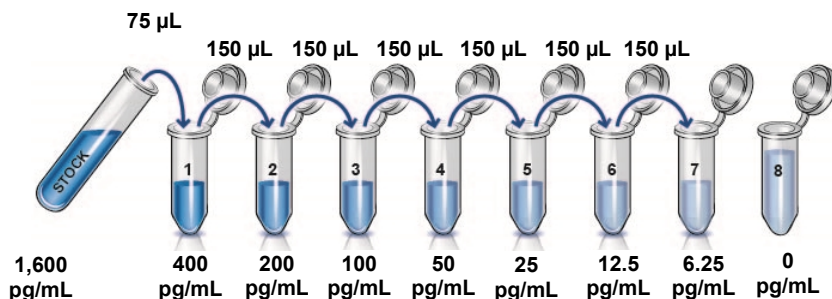
10.1 For measurements of **cell culture supernatant samples** follow this section to prepare the standard

10.1.1 Reconstitute the IL-17A standard by adding water by pipette.

10.1.2 Label eight tubes, Standards 1– 8.

10.1.3 Add 225 μL Sample Diluent NS into tube number 1 and 150 μL 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:

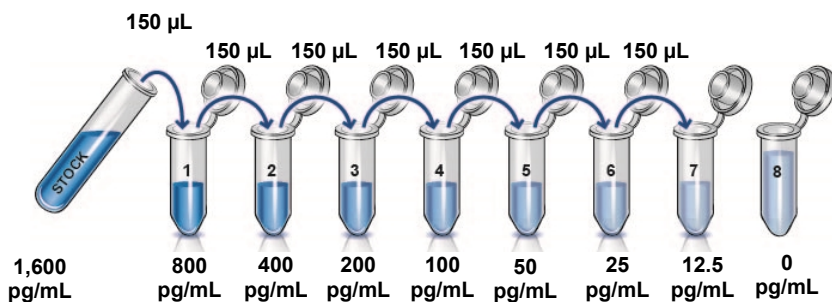


ASSAY PREPARATION

10.2 For measurements of **serum and plasma samples**, reconstitute the IL-17A standard by adding Sample Diluent 25BP by pipette.

10.2.1 Label eight tubes, Standards 1– 8 and 150 μ L Sample Diluent 25BP into each tube.

10.2.2 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
48 hours PHA+PMA treated EL4 cell culture supernatant	150 - 9,600X diluted
Mouse Plasma - Citrate	neat

11.1 Plasma

Collect plasma using citrate. Centrifuge samples at 2,000 x g for 10 minutes. Neat plasma samples can be assayed without dilution. If needed, dilute plasma samples into Sample Diluent 25BP 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. Neat serum samples can be assayed without dilution. If needed, dilute serum samples into Sample Diluent 25BP 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. Dilute supernatant samples at least 2X 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.

ASSAY PROCEDURE

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.

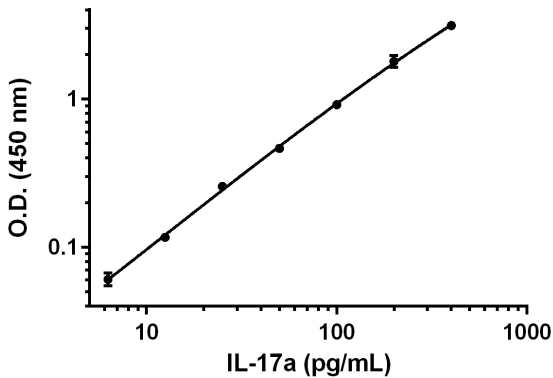
13.9 Analyze the data as described below.

14. CALCULATIONS

Subtract average zero standard from all readings. Average the duplicate readings of the positive control dilutions and plot against their concentrations. Draw the best smooth curve through these points to construct a standard curve. Most plate reader software 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 see which provides the most accurate (e.g. linear, semi-log, log/log, 4 parameter logistic). Interpolate protein concentrations for unknown samples from the standard curve plotted. Samples producing signals greater than that of the highest standard should be further diluted and reanalyzed, then multiplying the concentration found by the appropriate dilution factor.

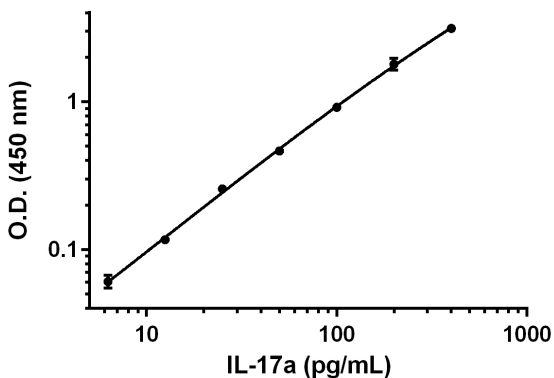
15. TYPICAL DATA

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	0.055	0.055	0.056
6.25	0.112	0.120	0.116
12.5	0.170	0.173	0.172
25	0.311	0.314	0.313
50	0.536	0.504	0.521
100	0.951	0.995	0.974
200	1.977	1.739	1.858
400	3.197	3.198	3.198

Figure 1. Example of IL-17A standard curve prepared in Sample Diluent NS. The IL-17A 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	0.054	0.051	0.053
12.5	0.127	0.128	0.128
25	0.203	0.201	0.203
50	0.376	0.363	0.370
100	0.705	0.708	0.707
200	1.376	1.388	1.383
400	2.710	2.717	2.714
800	3.804	3.774	3.790

Figure 2. Example of IL-17A standard curve prepared in Sample Diluent 25BP. The IL-17A 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	29	0.5 pg/mL
Sample Diluent 25BP	26	5.5 pg/mL

RECOVERY –

Three concentrations of mouse recombinant IL-17A 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% Mouse Serum	102.5	97.7 – 105.8
100% Mouse Plasma - Citrate	103.6	102.2 – 105.8
50% Cell Culture Media (with 10% Horse serum)	91.1	82.7 – 99.7

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 mouse IL-17A was measured in the following biological samples in a 2-fold dilution series. Sample dilutions were made in Sample Diluent NS.

Dilution Factor	Interpolated value	150X Diluted EL4 PHA+PMA-Stimulated Cell Culture Supernatant
Undiluted	pg/mL	426.42
	% Expected value	100
2	pg/mL	208.42
	% Expected value	97.8
4	pg/mL	97.71
	% Expected value	91.7
8	pg/mL	52.24
	% Expected value	98.0
16	pg/mL	28.53
	% Expected value	107
32	pg/mL	12.48
	% Expected value	93.7
64	pg/mL	7.64
	% Expected value	114.7

DATA ANALYSIS

Recombinant mouse IL-17A was spiked into the following biological samples and diluted in a 2-fold dilution series in Sample Diluent 25BP.

Dilution Factor	Interpolated value	100% Mouse Serum	100% Mouse Plasma (Citrate)
Undiluted	pg/mL	670.56	651.49
	% Expected value	100	100
2	pg/mL	423.38	322
	% Expected value	126.3	98.9
4	pg/mL	189.69	156.66
	% Expected value	113.2	96.2
8	pg/mL	98.20	88.08
	% Expected value	117.2	108.2
16	pg/mL	54.27	48.86
	% Expected value	129.5	120.0
32	pg/mL	27.18	22.38
	% Expected value	129.7	109.9

NL – Non-Linear

PRECISION –

Mean coefficient of variations of interpolated values from 3 concentrations of mouse recombinant IL17-a in neat mouse serum within the working range of the assay.

	Intra-Assay	Inter-Assay
n=	5	3
CV (%)	4.9	10.6

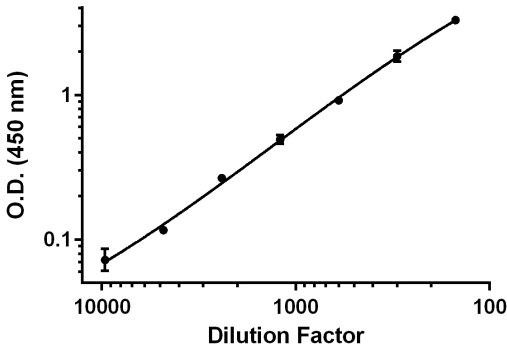


Figure 3. Demonstration of the linearity of dilution of the assay. Titration of PHA+PMA stimulated EL4 cell culture supernatant within the working range of the assay. EL4 cells were cultured in the presence of 1.5% phytohemagglutinin (PHA) and 10 ng/mL phorbol 12-myristate 13-acetate (PMA) for 48 hours. The cell culture supernatant was collected and measured in 2-fold dilution series. Background-subtracted data values (mean +/- SD, n = 2) are graphed.

Serum and Plasma. Neat pooled female mouse serum and plasma samples were measured for IL-17A concentrations using this kit. Means of interpolated sample values (n=5) are shown in pg of IL-17A per mL of serum/plasma in the table below. Note that IL-17A was not detectable (ND) in the neat pooled female mouse serum.

Sample Type	Mean (ng/mL)
Mouse serum	ND
Mouse Plasma - Citrate	13.0

17. ASSAY SPECIFICITY

This kit recognizes both native and recombinant mouse IL-17A protein in serum, plasma (citrate) and cell culture supernatant samples only.

Plasma-EDTA, plasma-heparin, milk, urine, saliva and cell and tissue extract samples have not been tested with this kit.

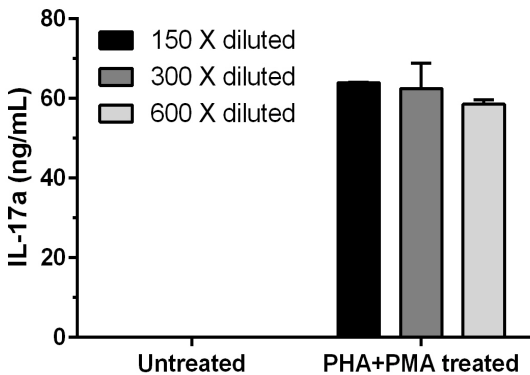


Figure 4. PHA+PMA treatment of EL4 cells stimulates secretion of IL-17A. EL4 cells were cultured in the absence or presence of 1.5% PHA and 10 ng/mL PMA for 48 hours. The cell culture supernatants were collected and their dilutions (as indicated) were measured with this kit. Interpolated concentrations of IL-17A adjusted for sample dilution are graphed in ng of IL-17A per mL of supernatant (mean \pm SD, n = 2). Note that IL-17A is not detectable in the untreated EL4 supernatant samples.

18. SPECIES REACTIVITY

This kit recognizes mouse and rat IL-17A proteins.

Rat species reactivity was determined only by measuring 2-fold dilutions of rat recombinant IL-17A protein (see below). Other species reactivity was not determined.

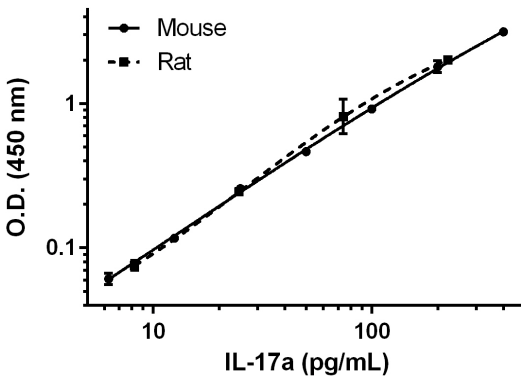


Figure 5. Titration of rat recombinant IL-17A within the working range of the assay. Two fold serial dilutions of mouse and rat recombinant IL-17A proteins were measured with this kit. Background-subtracted data values (mean +/- SD, n = 2) are graphed.

Please contact our Technical 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 © 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)