

Version 2 Last updated 19 November 2021

ab236715 Human IL-29 SimpleStep ELISA[®] Kit

For the quantitative measurement of IL-29 in human serum, plasma, and cell culture supernatants.

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	10
12. Plate Preparation	11
13. Assay Procedure	12
14. Calculations	14
15. Typical Data	15
16. Typical Sample Values	17
17. Assay Specificity	23
18. Species Reactivity	23
19. Troubleshooting	24
20. Notes	25
Technical Support	26

1. Overview

IL-29 *in vitro* SimpleStep ELISA® (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-29 protein in human serum, plasma, and cell culture supernatants.

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-29 (interleukin-29), also known as Interferon lambda-1, is a cytokine with antiviral, antitumor and immunomodulatory activities. It plays a critical role in the antiviral host defense, predominantly in the epithelial tissues. IL-29 is a ligand for the heterodimeric class II cytokine receptor composed of IL10RB and IFNLR1, and receptor engagement leads to the activation of the JAK/STAT signaling pathway resulting in the expression of IFN-stimulated genes, which mediate the antiviral state. IL-29 has a restricted receptor distribution and therefore restricted targets, it is primarily active in epithelial cells and this cell type-selective action is because of the epithelial cell-specific expression of its receptor IFNLR1. IL-29 exerts an immunomodulatory effect by up-regulating MHC class I antigen expression.

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
Human IL-29 Capture Antibody 10X	600 µL	+4°C
Human IL-29 Detector Antibody 10X	600 µL	+4°C
Human IL-29 Lyophilized Recombinant Protein	2 Vials	+4°C
Antibody Diluent 4BI	6 mL	+4°C
Wash Buffer PT 10X	20 mL	+4°C
TMB Development Solution	12 mL	+4°C
Stop Solution	12 mL	+4°C
Sample Diluent NS	50 mL	+4°C
Sample Diluent 50BP	20 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.

- 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.
- Sample Diluent BP may contain precipitate, this is normal. If precipitate is not dissolved by gentle mixing, the precipitate may be dissolved by gentle warming and mixing at 37°C for 10 minutes. If precipitate remains, gently spin down and avoid visible precipitates when pipetting.
- 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 10BP (for cell supernatants only):

Prepare **Sample Diluent 10BP** by diluting **Sample Diluent 50BP** with **Sample Diluent NS**. To make 10 mL Sample Diluent 10BP combine 2 mL Sample Diluent 50BP and 8 mL Sample Diluent NS. 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 4BI. 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 4BI. Mix thoroughly and gently.

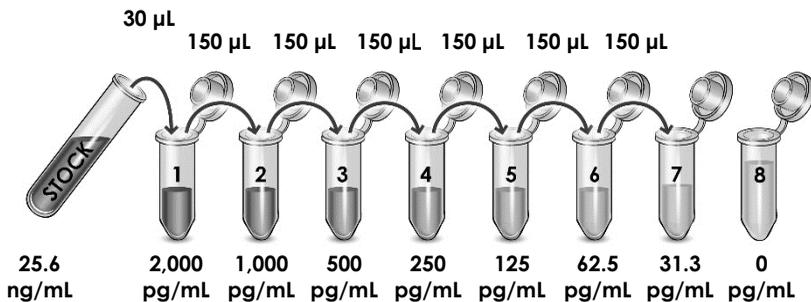
10. Standard Preparation

- 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).

IMPORTANT: If the protein standard vial has a volume identified on the label, reconstitute the IL-29 by adding that volume of Diluent indicated on the label. Alternatively, if the vial has a mass identified, reconstitute the IL-29 standard by adding 500 μ L Diluent.

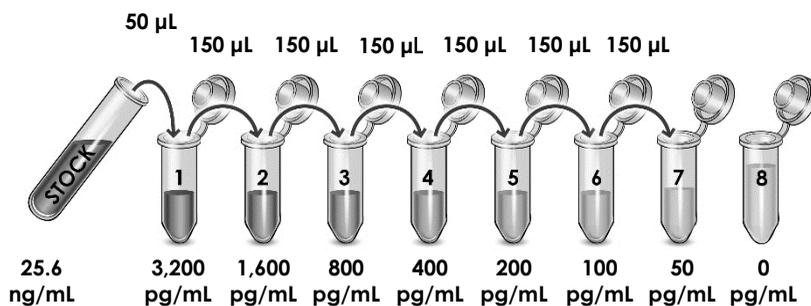
10.1 For cell culture supernatant samples follow these instructions:

- 10.1.1 Reconstitute the IL-29 standard sample by adding Sample Diluent 10BP. Hold at room temperature for 10 minutes and mix thoroughly and gently. This is the 25.6 ng/mL **Stock Standard Solution**.
- 10.1.2 Label eight tubes, Standards 1– 8.
- 10.1.3 Add 354 μ L of Sample Diluent 10BP into tube number 1 and 150 μ L of Sample Diluent 10BP 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:



10.2 For serum and plasma samples follow these instructions:

- 10.2.1 Reconstitute the IL-29 standard sample by adding Sample Diluent 50BP. Hold at room temperature for 10 minutes and mix thoroughly and gently. This is the 25.6 ng/mL **Stock Standard** Solution.
- 10.2.2 Label eight tubes, Standards 1– 8.
- 10.2.3 Add 350 μ L of Sample Diluent 50BP into tube number 1 and 150 μ L of Sample Diluent 50BP 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
Serum*	≤100%
Plasma – Citrate*	≤100%
Plasma – Heparin*	≤100%
Plasma – EDTA*	≤100%
IFN-alpha and Poly(I:C) Stimulated HUVEC Cell Culture Supernatant	12.5 – 50%

*Based on spiked sample

11.1 Plasma:

Collect plasma using citrate, EDTA or heparin. Centrifuge samples at 2,000 x g for 10 minutes. Samples may be assayed neat or dilute into Sample Diluent 50BP 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. Samples may be assayed neat or dilute into Sample Diluent 50BP 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 samples at least 2-fold into Sample Diluent 10BP and assay. Store un-diluted 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.
- 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 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. 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 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.

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.

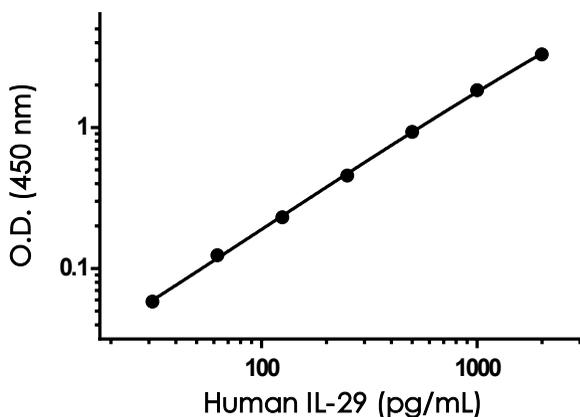
13.10 Analyze the data as described below.

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.

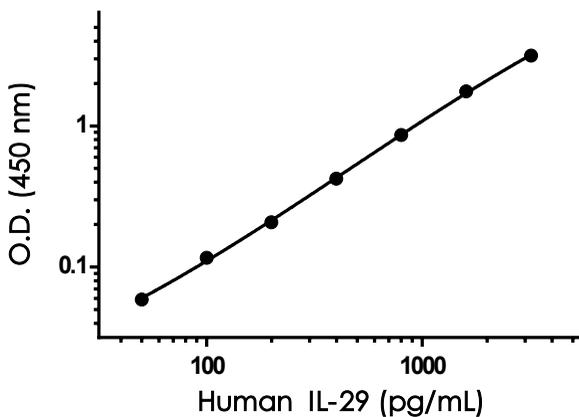
15. Typical Data

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	0.062	0.065	0.064
31.3	0.123	0.121	0.122
62.5	0.187	0.189	0.188
125	0.290	0.300	0.295
250	0.526	0.514	0.520
500	0.992	0.999	0.995
1,000	1.919	1.897	1.908
2,000	3.373	3.348	3.361

Figure 1. Example of human IL-29 standard curve in Sample Diluent 10BP. The IL-29 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	0.057	0.056	0.057
50	0.111	0.121	0.116
100	0.169	0.178	0.173
200	0.258	0.271	0.265
400	0.481	0.478	0.479
800	0.919	0.921	0.920
1,600	1.865	1.767	1.816
3,200	3.202	3.217	3.210

Figure 2. Example of human IL-29 standard curve in Sample Diluent 50BP. The IL-29 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 human IL-29. The NIBSC/WHO unclassified purified human IL-29 preparation 10/176 was evaluated in this kit.

The dose response curve of the unclassified standard IL-29 parallels the SimpleStep standard curve. To convert sample values obtained with the SimpleStep Human IL-29 kit to approximate NIBSC 10/176 units, use the equation below.

NIBSC (10/176) approximate value (IU/mL) = .03255 x SimpleStep Human IL-29 value (pg/mL).

17. Typical Sample Values

SENSITIVITY –

The MDD was determined by calculating the mean of zero standard replicates and adding 2 standard deviations then extrapolating the corresponding concentration.

Sample Diluent Buffer	n=	Minimal Detectable Dose
Sample Diluent 10BP	8	7.8 pg/mL
Sample Diluent 50BP	21	9.8 pg/mL

RECOVERY –

Three concentrations of IL-29 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 (%)
Neat Serum	111	107 – 115
Neat Plasma – Citrate	88	85 – 91
Neat Plasma – Heparin	117	116 – 117
Neat Plasma – EDTA	93	85 – 102
50% Unstimulated HUVEC Cell Culture Supernatant	99	95 – 102
50% Cell Culture Media*	106	102 – 109

*Media is RPMI 1640 containing 10% fetal calf 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.

Recombinant IL-29 was spiked into the following biological samples and diluted in a 2-fold dilution series. Serum and plasma samples are diluted in Sample Diluent 50BP. Cell culture media samples are diluted in Sample Diluent 10BP.

Dilution Factor	Interpolated value	100% Human Serum	100% Human Plasma (Citrate)	100% Human Plasma (Heparin)	100% Human Plasma (EDTA)	50% Cell Culture Media
Undiluted	pg/mL	1,986	1,520	1,819	1,840	1,593
	% Expected value	100	100	100	100	100
2	pg/mL	861.8	774.3	885.7	825.6	763.3
	% Expected value	87	102	97	90	96
4	pg/mL	417.3	385.3	406.6	416.9	372.8
	% Expected value	84	101	89	91	94
8	pg/mL	197.3	177.6	202.1	186.1	174.3
	% Expected value	90	93	89	81	88
16	pg/mL	108.6	80.07	91.39	NL	84.24
	% Expected value	88	84	80	NL	85

NL – Non-Linear

Neat pooled serum and plasma (EDTA, heparin, citrate) samples from healthy donors was measured in duplicate. All values were below the detectable range of the assay.

Neat serum from ten individual healthy human female/male donors was measured in duplicate. All values were below the detectable range of the assay.

PRECISION –

Mean coefficient of variations of interpolated values of IL-29 from three concentrations of IL-29 recombinant protein spiked into neat human serum within the working range of the assay.

	Intra- Assay	Inter- Assay
n =	5	3
CV (%)	2.8	2.5

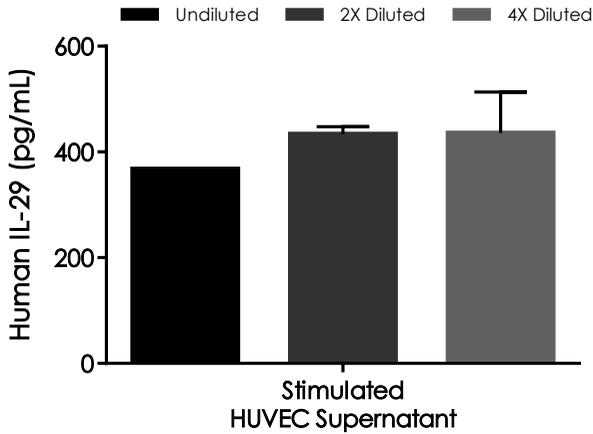


Figure 3. Interpolated concentrations of native IL-29 in IFN-alpha and Poly(I:C) stimulated HUVEC cell culture supernatant samples. The concentrations of IL-29 were measured in duplicates, interpolated from the IL-29 standard curves and corrected for sample dilution. Undiluted samples are as follows: 50% HUVEC supernatant. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean IL-29 concentration was determined to be 412.4 pg/mL in neat stimulated HUVEC cell culture supernatant.

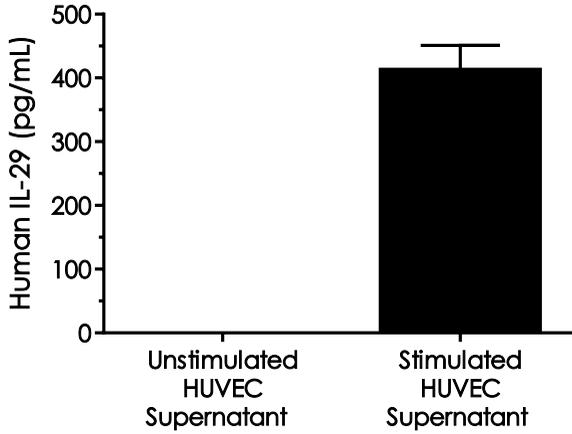


Figure 4. Comparison of IL-29 in unstimulated and IFN alpha and Poly(I:C) stimulated HUVEC cell culture supernatants. HUVEC cells were cultured for 20 hours in the presence or absence of 100 IU/mL human IFN alpha and 30 µg/mL Poly(I:C). The concentrations of IL-29 were measured in three different dilutions of the supernatant samples in duplicates and interpolated from the IL-29 standard curve. The interpolated values are plotted (mean +/- SD, n=3). The mean IL-29 concentration was determined to be 412.4 pg/mL in stimulated HUVEC cell culture supernatant, and undetectable in unstimulated HUVEC cell culture supernatant and RPMI1640 media containing 10% FBS (not shown).

18. Assay Specificity

This kit recognizes both native and recombinant human IL-29 protein in serum, plasma, and cell culture supernatant samples only.

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

CROSS REACTIVITY

Recombinant human IL-28A was prepared at 1 ng/mL and assayed for cross reactivity. No cross-reactivity was observed.

INTERFERENCE

Recombinant human IL-28A was prepared at 50 ng/mL and 1 ng/mL and tested for interference with 1 ng/mL recombinant human IL-29. No interference with was observed.

19. Species Reactivity

This kit recognizes human IL-29 protein.

Other species reactivity not determined.

Please contact our Technical Support team for more information.

20. Troubleshooting

Problem	Reason	Solution
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.

21. 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)