



**ab183364 –  
Resistin (RETN) Human  
SimpleStep ELISA<sup>®</sup> Kit**

Instructions for Use

For the quantitative measurement of Resistin (RETN) 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

---

## INTRODUCTION

- 1. BACKGROUND 2
- 2. ASSAY SUMMARY 4

## GENERAL INFORMATION

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

## ASSAY PREPARATION

- 9. REAGENT PREPARATION 8
- 10. STANDARD PREPARATION 9
- 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 20
- 18. SPECIES REACTIVITY 20

## RESOURCES

- 19. TROUBLESHOOTING 21
- 20. NOTES 22

## 1. BACKGROUND

Abcam's Resistin (RETN) *in vitro* SimpleStep ELISA® (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Resistin 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 substrate 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.

Resistin, otherwise known as Found in Inflammatory Zone 3 (FIZZ3), and Adipocyte Secreted Factor (ADSF), is a 114 amino-acid cysteine-rich protein. It can be found as a high molecular weight hexamer or as a more active, lower molecular weight complex. Resistin is the best known member of the family of Resistin-like Molecules (RELMs), which include RELM-alpha/FIZZ1 and RELM-beta/FIZZE 2, both described only in rodents. Furthermore, Resistin has been categorized as part of the group of adipokines hormones, which are known to be involved in the development of obesity, insulin resistance, diabetes, inflammation, autoimmune disease and metabolic syndrome.

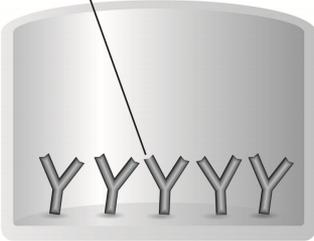
In Humans, Resistin is a pro-inflammatory molecule primarily secreted by macrophages and mononuclear bone marrow cells in response to

inflammation, hyperglycemia, and growth hormones. IL-1, IL-6, TNF-alpha as well as lipopolysaccharide (LPS) are strong inducers of resistin as observed in animal models of hepatic inflammation and necrosis. High levels of Resistin are also observed in patients with pancreatitis and with systemic lupus erythematosus.

Beyond its role as a pro-inflammatory molecule, Resistin has been found to have a metabolic function due to its expression in other tissues such as white adipose tissue, placenta, uterus, skeletal muscle, GI tract, spleen, thyroid gland and pancreas. Metabolically, elevated levels of Resistin are associated with glucose intolerance, hyperinsulinemia, and insulin resistance. These effects are believed to be consequential to dephosphorylation of AMPK in the liver and overexpression of Suppressor of cytokine signaling 3 (SOCS-3) in the pancreatic beta cells. Furthermore, the levels of Resistin in serum have been found to positively correlate with increased body fat content, which may be explained by its ability to induce adipocyte proliferation and angiogenesis. Despite the effects of Resistin on insulin resistance, glycemic levels and body fat content, the role of Resistin linking obesity to diabetes continues to be controversial.

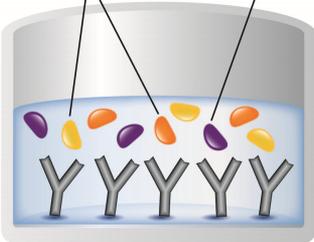
## 2. ASSAY SUMMARY

Immobilization Antibody



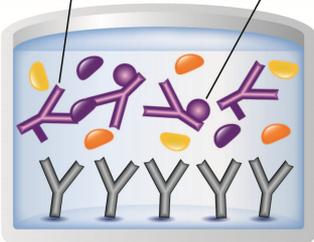
Remove appropriate number of antibody coated well strips. Equilibrate all reagents to room temperature. Prepare all reagents, samples, and standards as instructed.

Matrix Proteins Target Analyte



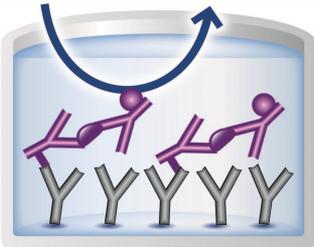
Add standard or sample to appropriate wells.

Capture Antibody Detector Antibody



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

Substrate Color Development



Aspirate and wash each well. Add TMB Substrate 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 sections 9 & 10.

### 5. MATERIALS SUPPLIED

Item	Amount	Storage Condition (Before Preparation)
10X Resistin Capture Antibody	600 µL	+2-8°C
10X Resistin Detector Antibody	600 µL	+2-8°C
Resistin Human 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 Substrate	12 mL	+2-8°C
Stop Solution	12 mL	+2-8°C
Sample Diluent NS	50 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<sub>2</sub>PO<sub>4</sub>, 8 mM Na<sub>2</sub>HPO<sub>4</sub>, 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.

## GENERAL INFORMATION

- 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 5X Cell Extraction Buffer contains phosphatase inhibitors. Protease inhibitors can be added if required.
- The provided 50X Cell Extraction Enhancer Solution may precipitate when stored at +4°C. To dissolve, warm briefly at +37°C and mix gently. The 50X Cell Extraction Enhancer Solution can be stored at room temperature to avoid precipitation.
- **To avoid high background always add samples or standards to the well before the addition of the antibody cocktail.**
- **This kit is sold based on number of tests. A 'test' simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.**

## 9. REAGENT PREPARATION

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

### 9.1 **1X 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**

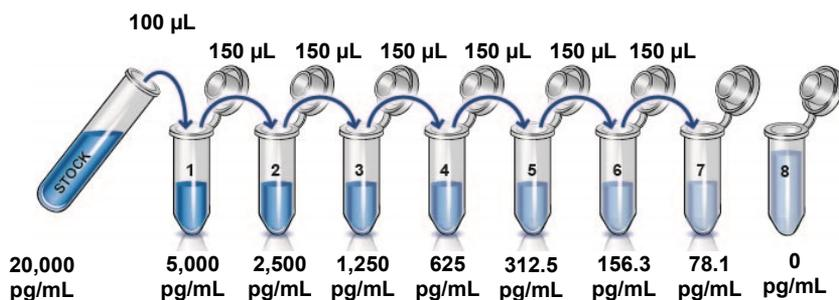
Prepare Antibody Cocktail by diluting the capture and detector antibodies in Antibody Diluent CPI2. To make 3 mL of the Antibody Cocktail combine 300  $\mu$ L 10X Capture Antibody and 300  $\mu$ L 10X Detector Antibody with 2.4 mL Antibody Diluent CPI2. Mix thoroughly and gently.

## 10. STANDARD PREPARATION

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

The following 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 Resistin standard by adding that volume of Sample Diluent NS indicated on the label. Alternatively, if the vial has a mass identified, reconstitute the Resistin standard by adding 500  $\mu\text{L}$  Sample Diluent NS. Hold at room temperature for 10 minutes and mix thoroughly and gently. This is the 20,000  $\text{pg/mL}$  **Stock Standard Solution** (see table below).
- 10.2 Label eight tubes with numbers 1 – 8.
- 10.3 Add 300  $\mu\text{L}$  Sample Diluent NS into tube #1 and 150  $\mu\text{L}$  Sample Diluent NS into tube numbers 2-8.
- 10.4 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 (%)
Normal Human Serum	12.5 – 100
Normal Human Plasma - Citrate	12.5 – 100
Normal Human Plasma - Heparin	12.5 – 100
Normal Human Plasma - EDTA	12.5 – 100
Cell culture media	12.5 – 100

### 11.1 Plasma

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

### 11.2 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. Dilute samples into Sample Diluent NS 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. If necessary, dilute supernatants in non-reactive cell culture media 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 Substrate 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  $\mu$ L Stop Solution to each well and recording the OD at 450 nm.*

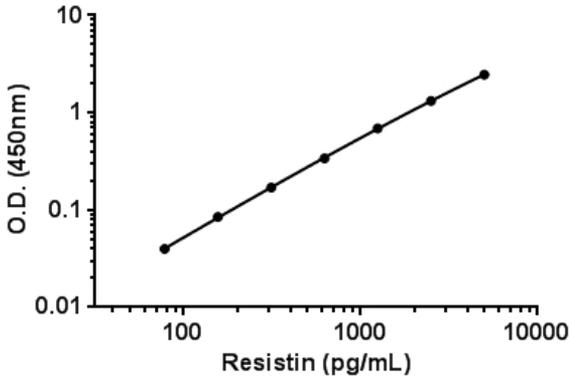
13.9 Analyze the data as described below.

## 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.  
  
*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.

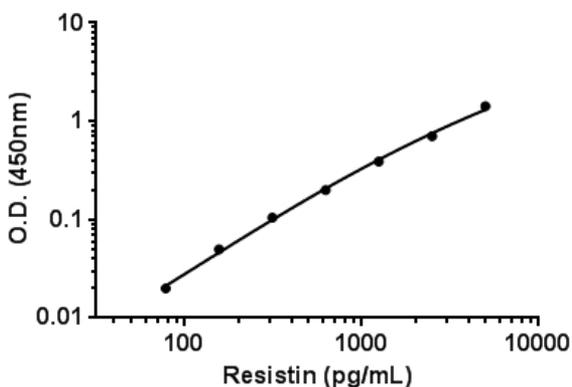
## 15. TYPICAL DATA

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



Standard Curve Measurements			
Conc. (pg/mL)	O.D. 450 nm		Mean O.D.
	1	2	
0	0.081	0.081	0.081
78.1	0.116	0.116	0.116
156.3	0.167	0.162	0.165
312.5	0.248	0.248	0.248
625	0.423	0.418	0.420
1,250	0.766	0.765	0.765
2,500	1.391	1.422	1.406
5,000	2.542	2.545	2.543

**Figure 1.** Example of Resistin standard curve in Sample Diluent NS. The Resistin 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	0.052	0.052	0.052
78.1	0.079	0.079	0.079
156.3	0.105	0.109	0.107
312.5	0.165	0.164	0.165
625	0.260	0.258	0.259
1,250	0.446	0.454	0.450
2,500	0.788	0.740	0.764
5,000	1.478	1.492	1.485

**Figure 2.** Example of Resistin standard curve in cell culture media. The Resistin 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 (MDD) dose is 24 pg/mL. The MDD was determined by calculating the mean of zero standard replicates (n=40) and adding 2 standard deviations then extrapolating the corresponding concentrations.

### **RECOVERY –**

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

<b>Sample Type</b>	<b>Average % Recovery</b>	<b>Range (%)</b>
100% Normal Human Serum	94	88 - 100
100% Normal Human Plasma - Citrate	78	68 - 90
100% Normal Human Plasma - EDTA	96	88 - 104
100% Normal Human Plasma - Heparin	95	84 - 110
Cell Culture Media	106	106 - 107

## 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.

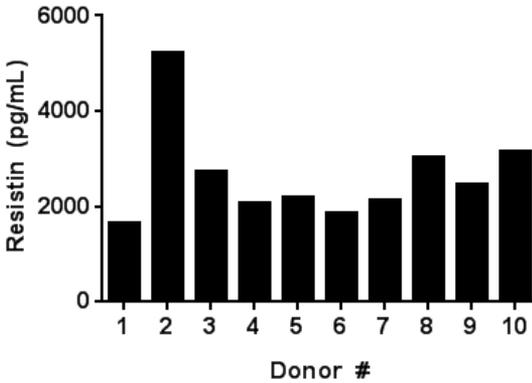
Recombinant Resistin was spiked into the following biological samples and diluted in a 2-fold dilution series in Sample Diluent NS.

Dilution Factor	Interpolated value	100% Human Serum	100% Human Plasma (Citrate)	100% Human Plasma (EDTA)	100% Human Plasma (Heparin)	100% cell culture media
Undiluted	pg/mL	2,567.8	2,421.8	2,688.0	2,789.9	3161.4
	<b>% Expected value</b>	<b>100</b>	<b>100</b>	<b>100</b>	<b>100</b>	<b>100</b>
2	pg/mL	1,302.8	1,387.9	1,435.2	1,387.9	1,499.4
	<b>% Expected value</b>	<b>101</b>	<b>115</b>	<b>107</b>	<b>108</b>	<b>95</b>
4	pg/mL	701.8	584.0	674.5	583.9	630.3
	<b>% Expected value</b>	<b>109</b>	<b>96</b>	<b>100</b>	<b>120</b>	<b>80</b>
8	pg/mL	362.5	337.5	375.3	442.4	365.6
	<b>% Expected value</b>	<b>113</b>	<b>111</b>	<b>112</b>	<b>142</b>	<b>93</b>

## PRECISION –

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

	Intra-Assay	Inter-Assay
n=	9	3
CV (%)	3	6



**Figure 3.** Levels of Resistin in normal serum. The levels of native Resistin in serum samples were tested from ten individual healthy donors. Levels were interpolated from the standard curve in Sample Diluent NS. The mean level of Resistin was 2.6 ng/mL with a range of 1.6 to 5.3 ng/mL and a standard deviation of 1 ng/mL.

### **17. ASSAY SPECIFICITY**

This assay recognizes native and recombinant human Resistin protein.

Cross-reactivity was tested with human Leptin recombinant protein prepared at 50ng/mL in Sample Diluent NS. No significant cross-reactivity was observed with OD deviation from background of 0.019. Similarly, recombinant human Resistin protein was assayed at 1,000 pg/mL in the presence and absence of 50 ng/mL of human Leptin to determine interference. Recovery of Resistin was observed at 103% with a standard deviation of 1%.

### **18. SPECIES REACTIVITY**

This kit recognizes both native and recombinant human Resistin protein in the serum, plasma, and cell culture supernatants only. Mouse and rat serum and plasma samples have not been tested with this kit.

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 substrate solution protected from light.

20. NOTES

## **Technical Support**

Copyright © 2024 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](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)