# Sigma-Aldrich

User Guide

# Mouse IL-17B ELISA Kit

For Serum, Plasma, Cell Culture Supernatants

#### **RAB0264**

#### Introduction

The Mouse IL-17B ELISA kit is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of mouse IL-17B in serum (mouse IL-17B concentration is low in normal serum/plasma and may not be detectable in this assay), plasma and cell culture supernatants. This assay employs an antibody specific for mouse IL-17B coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-17B present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse IL-17B antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of IL-17B bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

## Storage

The entire kit may be stored at -20°C for up to 1 year from date of shipment. Avoid repeated freeze thaw cycles. The kit may be stored at 4°C for up to 6 months. For extended storage, it is recommended to store at -80°C.

## Components

- Mouse IL-17B Microplate RAB0264A-EA: 96 wells (12 strips x 8 wells) coated with anti-Mouse IL-17B. Stable in storage 1 month at 4 °C once opened. Return unused wells to the pouch containing desiccant pack, reseal along the edge.
- Mouse IL-17B Standard Protein RAB0264C-1VL: 2 vials of Mouse IL-17B. 1 vial is enough to run each standard in duplicate. Stable in storage 1 week at -80 °C once opened.
- Mouse IL-17B Detection Antibody RAB0264D-1VL: 2 vials of biotinylated anti-Mouse IL-17B. Each vial is enough to assay half the microplate. Stable in storage 5 days at 4 °C once opened.
- Wash Buffer RABWASH4: 25 mL of 20X concentrated solution. Stable in storage 1 month at 4 °C once opened.
- HRP-Streptavidin RABHRP5: 200 μL 200X concentrated HRP-conjugated streptavidin. Do not store and reuse.
- TMB One-Step Substrate Reagent RABTMB3: 12 mL of 3,3',5,5'-tetramethylbenzidine (TMB) in buffer solution.
- Stop Solution RABSTOP3: 8 mL of 0.2 M sulfuric acid.
- 5X Assay Diluent D RABELADD-15ML: 15 mL of 5X concentrated buffer. Stable in storage 1 month at 4°C once opened.
- 5X Assay Diluent B RABELADE-15ML: 15 mL of 5X concentrated buffer. Stable in storage 1 month at 4°C once opened.



## Additional Materials Required (But Not Provided)

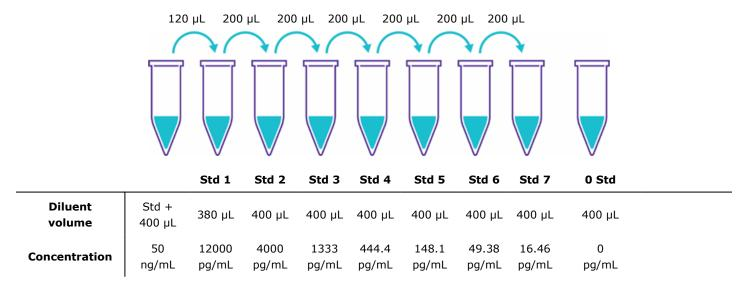
- Microplate reader capable of measuring absorbance at 450 nm.
- Precision pipettes to deliver 2 µl to 1 mL volumes.
- Adjustable 1-25 mL pipettes for reagent preparation.
- 100 mL and I liter graduated cylinders.
- · Absorbent paper.
- Distilled or deionized water.
- Log-log graph paper or computer and software for ELISA data analysis.
- Tubes to prepare standard or sample dilutions.

## Reagent Preparation

- 1. Bring all reagents and samples to room temperature (18-25 °C) before use.
- 2. 5X Assay Diluent D and 5X Assay Diluent B should be diluted 5-fold with deionized or distilled water before use.
- 3. Sample dilution: 1X Assay Diluent D should be used for dilution of serum, plasma, and cell culture supernatant samples. The suggested dilution for normal serum/plasma is 2-fold.

**Note:** Levels of IL-17B may vary between different samples. Optimal dilution factors for each sample must be determined by the investigator.

4. Preparation of standard: Briefly spin a vial of Standard Protein. Add 400 μL 1X Assay Diluent D into the Standard Protein vial to prepare a 50 ng/mL standard solution. Dissolve the powder thoroughly by a gentle mix. Add 120 μL IL-17B standard from the vial of Standard Protein, into a tube with 380 μL 1X Assay Diluent D to prepare a 12,000 pg/mL standard solution. Pipette 400 μL 1X Assay Diluent D into each tube. Use the stock standard solution to produce a dilution series (shown below). Mix each tube thoroughly before the next transfer. 1X Assay Diluent D serves as the zero standard (0 pg/mL).



- 5. If the Wash Buffer (20X) contains visible crystals, warm to room temperature and mix gently until dissolved. Dilute 20 mL of Wash Buffer Concentrate into deionized or distilled water to yield 400 mL of 1X Wash Buffer.
- 6. Briefly spin the Detection Antibody vial before use. Add 100 μL of 1X Assay Diluent B into the vial to prepare a detection antibody concentrate. Pipette up and down to mix gently (the concentrate can be stored at 4 °C for 5 days). The detection antibody concentrate should be diluted 80-fold with 1X Assay Diluent B and used in step 5 of the assay procedure.
- 7. Briefly spin the HRP-Streptavidin concentrate vial and pipette up and down to mix gently before use, as precipitates may form during storage. HRP-Streptavidin concentrate should be diluted 200-fold with 1X Assay Diluent B.

For example: Briefly spin the HRP-Streptavidin vial and pipette up and down to mix gently. Add 50  $\mu$ L of HRP-Streptavidin concentrate into a tube with 10 mL 1X Assay Diluent B to prepare a 200-fold diluted HRP-Streptavidin solution (don't store the diluted solution for next day use). Mix well.

## **Assay Procedure**

- 1. Bring all reagents and samples to room temperature (18-25 °C) before use. It is recommended that all standards and samples be run at least in duplicate.
- 2. Label removable 8-well strips as appropriate for your experiment.
- 3. Add 100  $\mu$ L of each standard (see Reagent Preparation step 4) and sample into appropriate wells. Cover wells and incubate for 2.5 hours at room temperature with gentle shaking.
- 4. Discard the solution and wash 4 times with 1X Wash Solution. Wash by filling each well with Wash Buffer (300  $\mu$ L) using a multi-channel Pipette or autowasher. Complete removal of liquid at each step is essential to good performance. After the last wash, remove any remaining Wash Buffer by aspirating or decanting. Invert the plate and blot it against clean paper towels.
- 5. Add 100  $\mu$ L of 1X prepared biotinylated antibody (Reagent Preparation step 6) to each well. Incubate for 1 hour at room temperature with gentle shaking.
- 6. Discard the solution. Repeat the wash as in step 4.
- 7. Add 100  $\mu$ L of prepared Streptavidin solution (see Reagent Preparation step 7) to each well. Incubate for 45 minutes at room temperature with gentle shaking.
- 8. Discard the solution. Repeat the wash as in step 4.
- 9. Add 100  $\mu$ L of TMB One-Step Substrate Reagent to each well. Incubate for 30 minutes at room temperature in the dark with gentle shaking.
- 10. Add 50 µL of Stop Solution to each well. Read at 450 nm immediately.

## Assay Procedure Summary

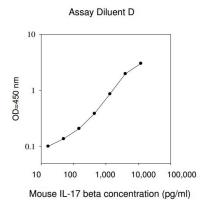
- 1. Prepare all reagents, samples and standards as instructed.
- 2. Add 100 µL standard or sample to each well. Incubate 2.5 hours at room temperature.
- 3. Add 100 µL prepared biotin antibody to each well. Incubate 1 hour at room temperature.
- 4. Add 100 μL prepared Streptavidin solution. Incubate 45 minutes at room temperature.
- 5. Add 100 μL TMB One-Step Substrate Reagent to each well. Incubate 30 minutes at room temperature.
- 6. Add 50 µL Stop Solution to each well. Read at 450 nm immediately.

#### Calculation of Results

Calculate the mean absorbance for each set of duplicate standards, controls and samples, and subtract the average zero standard optical density. Plot the standard curve on log-log graph paper or using Sigma plot software, with standard concentration on the x-axis and absorbance on the y-axis. Draw the best-fit straight line through the standard points.

#### Typical Data

These standard curves are for demonstration only. A standard curve must be run with each assay.



#### Sensitivity

The minimum detectable dose of Mouse IL-17B was determined to be 12 pg/mL.

Minimum detectable dose is defined as the analyte concentration resulting in an absorbance that is 2 standard deviations higher than that of the blank (diluent buffer).

## Spiking & Recovery

Recovery was determined by spiking various levels of Mouse IL-17B into the sample types listed below. Mean recoveries are as follows:

Sample Type	Average % Recovery	Range (%)
Serum	119	104-128
Plasma	129	92-155
Cell Culture Supernatants	149	139-161

#### Linearity

Sample Type		Serum	Plasma	Cell Culture Supernatants
1:2	Average % of Expected	109	108	106
1.2	Range (%)	95-117	98-117	94-113
1:4	Average % of Expected Range	105	113	106
1.4	Range (%)	94-115	102-124	94-104

#### Reproducibility

Intra-Assay CV%: <10% Inter-Assay CV%: <12%

## Specificity

This ELISA kit shows no cross-reactivity with the following cytokines tested: Mouse CD30, L CD30, T CD40, CRG-2, CTACK, CXCL16, Eotaxin , Eotaxin-2, Fas Ligand, Fractalkine, GCSF, IFN- gamma, IGFBP-3, IGFBP-5, IGFBP-6, IL-1 alpha, IL-1 beta, IL-13, KC, Leptin R, LEPTIN(OB), LIX, L-Selectin, Lymphotactin, MCP-5, M-CSF, MIG, MIP- 1alpha, MIP-1 gamma, MIP-2, MIP-3 beta, MIP-3 alpha, PF-4, P-Selectin, SCF, SDF-1 alpha, TARC, TCA-3, TECK, TIMP-1, TNF RI, TPO, VEGF.

## Troubleshooting Guide

Problem	Cause	Solution		
	Inaccurate pipetting	Check pipettes		
Poor standard curve	Improper standard dilution	Briefly centrifuge the standard protein and dissolve the powder thoroughly by gently mixing		
Low signal	Improper preparation of standard and/or biotinylated antibody	Briefly spin down vials before opening. Dissolve the powder thoroughly.		
	Too brief incubation times	Ensure sufficient incubation time. Assay procedure step 3 may be done overnight at 4 °C with gentle shaking.		
		Note: may increase overall signals including background.		
	Inadequate reagent volumes or improper dilution	Check pipettes and ensure correct preparation		
Large CV	Inaccurate pipetting	Check pipettes		
	Air bubblies in wells	Remove bubbles in wells		
High background	Plate is insufficiently washed	Review the manual for proper wash. If using a plate washer, ensure that all ports are unobstructed.		
	Contaminated wash buffer	Make fresh wash buffer		
Low sensitivity	Improper storage of the ELISA kit	Store your standard at -70 °C after reconstitution, others at 4 °C. Keep substrate solution protected from light.		
	Stop Solution	Add stop solution to each well before reading plate		

#### **Notice**

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

#### **Technical Assistance**

Visit the tech service page at SigmaAldrich.com/techservice.

#### Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at SigmaAldrich.com/terms.

#### **Contact Information**

For the location of the office nearest you, go to SigmaAldrich.com/offices.



RAB0264pis Rev 08/24