

### Mouse IL-17B ELISA Kit ab100703

#### 画像数 1

#### 製品の概要

製品名	Mouse IL-17B ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 12 pg/ml
検出範囲	16.46 pg/ml - 12000 pg/ml
添加回収試験	> 100 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	149.4	139% - 161%
Serum	118.9	104% - 128%
Plasma	129.3	92% - 155%

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

製品の概要 Abcam's IL-17B Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse IL-17B in serum (mouse IL-17B concentration is pretty low in normal serum/plasma, it may not be detected 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.

試験プラットフォーム Microplate

## 製品の特性

### 保存方法

Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
Biotinylated anti-mouse IL-17B	2 vials
IL-17B Microplate (12 x 8 wells)	1 unit
Recombinant Mouse IL-17B Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

### 機能

Stimulates the release of tumor necrosis factor alpha and IL-1-beta from the monocytic cell line THP-1.

### 組織特異性

Expressed in adult pancreas, small intestine, stomach, spinal cord and testis. Less pronounced expression in prostate, colon mucosal lining, and ovary.

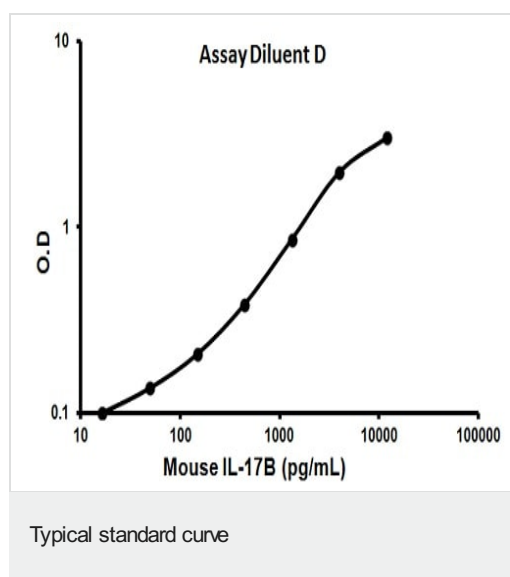
### 配列類似性

Belongs to the IL-17 family.

### 細胞内局在

Secreted.

## 画像



Representative standard curve using ab100703

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

## **Our Abpromise to you: Quality guaranteed and expert technical support**

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
  
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.co.jp/abpromise> or contact our technical team.

## **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors