

Mouse IL-17B ELISA Kit ab100703

1 图像

概述

| | |
|------|---|
| 产品名称 | 小鼠IL-17B ELISA试剂盒 |
| 检测方法 | Colorimetric |
| 样品类型 | Cell culture supernatant, Serum, Plasma |
| 检测类型 | Sandwich (quantitative) |
| 灵敏度 | < 12 pg/ml |
| 范围 | 16.46 pg/ml - 12000 pg/ml |
| 回收率 | > 100 % |

特定样本回收率

| 样品类型 | 平均 % | 范围 |
|--------------------------|-------|-------------|
| Cell culture supernatant | 149.4 | 139% - 161% |
| Serum | 118.9 | 104% - 128% |
| Plasma | 129.3 | 92% - 155% |

实验步骤 Multiple steps standard assay

种属反应性 与反应: Mouse

产品概述 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.

平台 Microplate

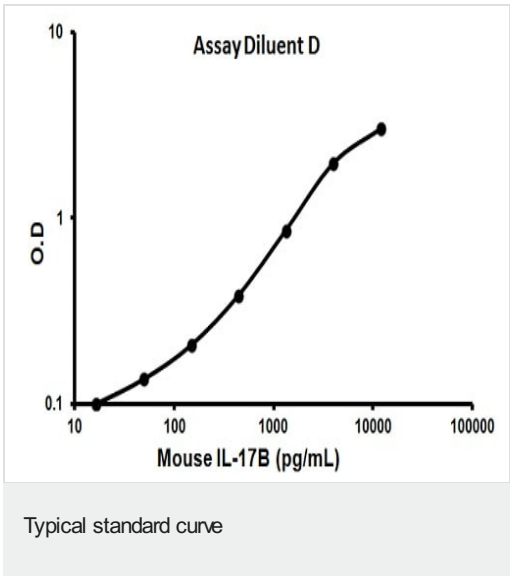
性能

存放说明 Store at -20°C. Please refer to protocols.

| 组件 | 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 |

| | |
|-------|---|
| 功能 | Stimulates the release of tumor necrosis factor alpha and IL-1-beta from the monocytic cell line THP-1. |
| 组织特异性 | Expressed in adult pancreas, small intestine, stomach, spinal cord and testis. Less pronounced expression in prostate, colon mucosal lining, and ovary. |
| 序列相似性 | Belongs to the IL-17 family. |
| 细胞定位 | Secreted. |

图片



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.cn/abpromise> or contact our technical team.

Terms and conditions

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