

16 References 画像数 2

製品の概要

製品名	Mouse IL-17 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	< 6 pg/ml
検出範囲	6.1 pg/ml - 600 pg/ml
添加回収試験	99 %

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	99.62	85% - 105%
Serum	101.38	87% - 107%
Plasma	97.47	84% - 103%

ステップ	Multiple steps standard assay
種交差性	交差種: Mouse
製品の概要	<p>Abcam's IL-17 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of mouse IL-17 in serum, plasma and cell culture supernatants.</p> <p>This assay employs an antibody specific for mouse IL-17 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-17 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse IL-17 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-17 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p>
試験プラットフォーム	Microplate

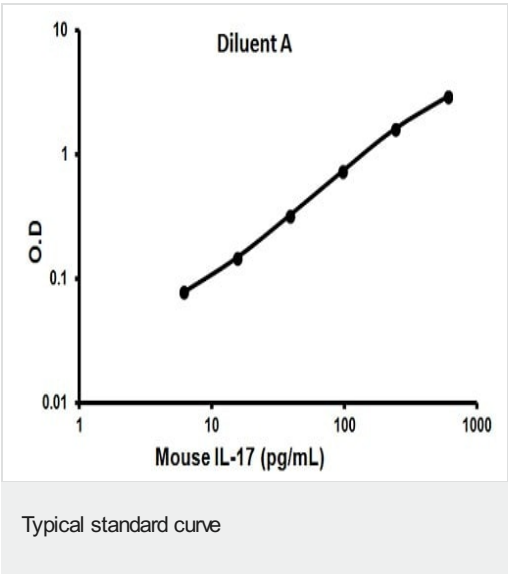
製品の特 性

保存方法Store at -20°C. Please refer to protocols.

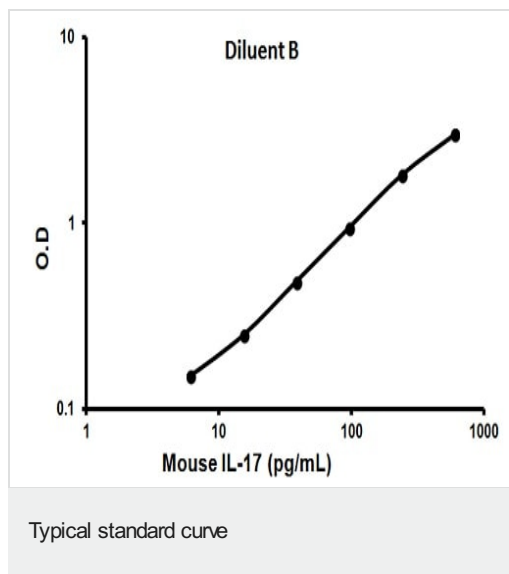
内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
700X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Mouse IL-17	2 vials
IL-17 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse IL-17 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

機能	Induces stromal cells to produce proinflammatory and hematopoietic cytokines. Enhances the surface expression of the intracellular adhesion molecule-1 (ICAM-1) in fibroblasts.
組織特異性	Restricted to activated memory T-cells.
配列類似性	Belongs to the IL-17 family.
翻訳後修飾	Found both in glycosylated and nonglycosylated forms.
細胞内局在	Secreted.

画像



Representative standard curve using ab100702



Representative standard curve using ab100702

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