

Mouse IL-21 ELISA Kit ab100707

画像数 2

製品の概要

| | |
|---------|----------------------------------|
| 製品名 | Mouse IL-21 ELISA Kit |
| 検出方法 | Colorimetric |
| サンプルの種類 | Cell culture supernatant, Plasma |
| アッセイタイプ | Sandwich (quantitative) |
| 検出感度 | < 6 pg/ml |
| 検出範囲 | 8.19 pg/ml - 2000 pg/ml |
| 添加回収試験 | 89 % |

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 | 98.65 | 87% - 109% |
| Plasma | 77.38 | 68% - 87% |

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

製品の概要

Abcam's IL-21 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse IL-21 in plasma and cell culture supernatants. (Mouse IL-21 concentration is pretty low in normal plasma, it may not be detected in this assay).

This assay employs an antibody specific for mouse IL-21 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-21 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse IL-21 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-21 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

We have not been able to detect endogenous Mouse IL-21 in normal serum with ab100707, only in serum spiked with Mouse IL-21.

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 |
| Assay Diluent A | 1 x 30ml |
| Biotinylated anti-mouse IL-21 | 2 vials |
| IL-21 Microplate (12 x 8 wells) | 1 unit |
| Recombinant mouse IL-21 Standard (lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

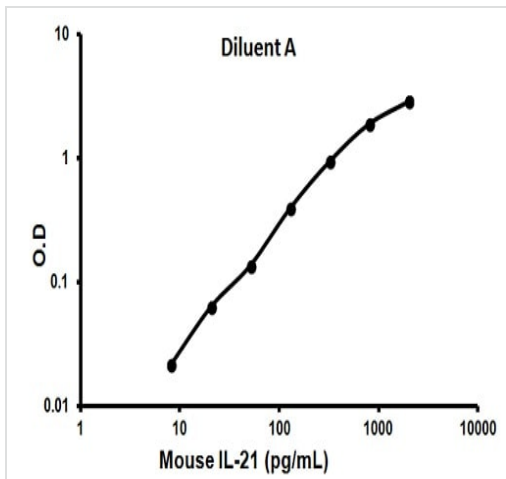
関連性

A novel cytokine related to IL-2 and IL-15 was recently identified and designated IL-21. IL-21 has been found to be a powerful growth factor for naive B cells. The receptor for IL-21 (IL-21R, also termed NILR for novel Interleukin receptor) is a new member of the class I cytokine receptor family. IL-21R forms a complex with the common cytokine receptor γ chain, γc , and mediates IL-21 signaling. Both IL-21R and the γc are necessary for the IL-21 function. IL-21 and its receptor activate JAK/STAT signaling pathway. IL-21R is expressed in spleen, thymus, natural killer (NK), T and B cell lines. IL-21 plays a role in the proliferation and maturation of NK, B and T cell populations.

細胞内局在

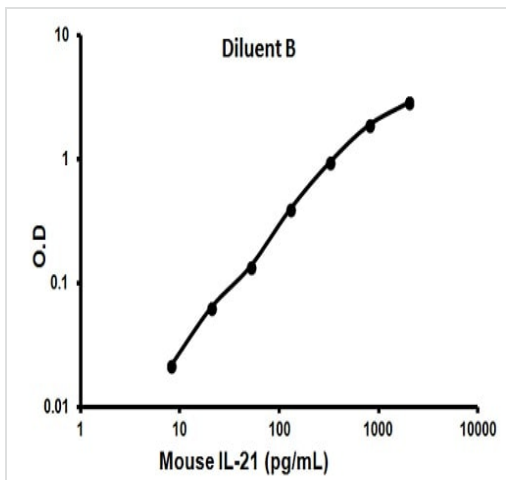
Secreted

画像



Representative standard curve using ab100707

Typical standard curve



Representative standard curve using ab100707

Typical standard curve

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