

Mouse IL-21 ELISA Kit ab100707

2 图像

概述

产品名称	小鼠IL-21 ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 6 pg/ml
范围	8.19 pg/ml - 2000 pg/ml
回收率	89 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	98.65	87% - 109%
Plasma	77.38	68% - 87%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Mouse

产品概述 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.

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
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

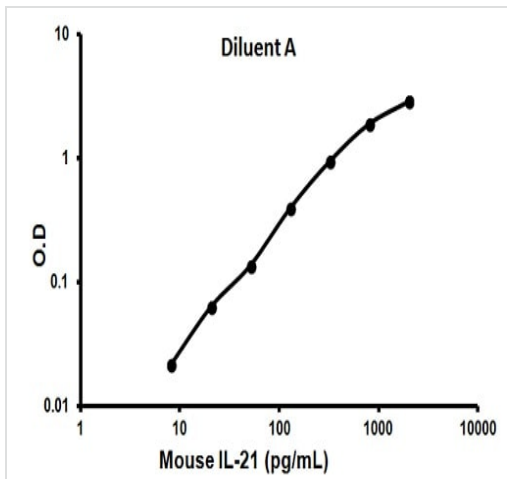
相关性

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.

细胞定位

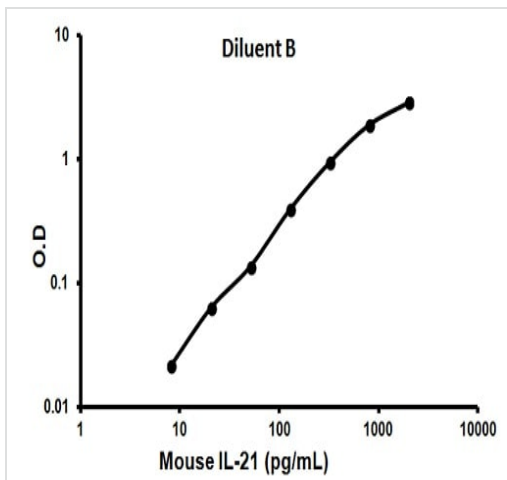
Secreted

图片



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.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