

Human IL-18 R beta ELISA Kit ab113325

1 图像

概述

产品名称	人IL-18 R beta 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	12.3 pg/ml - 3000 pg/ml
回收率	77 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	75.78	68% - 85%
Serum	80.9	71% - 89%
Plasma	74.98	67% - 83%

实验步骤 Multiple steps standard assay

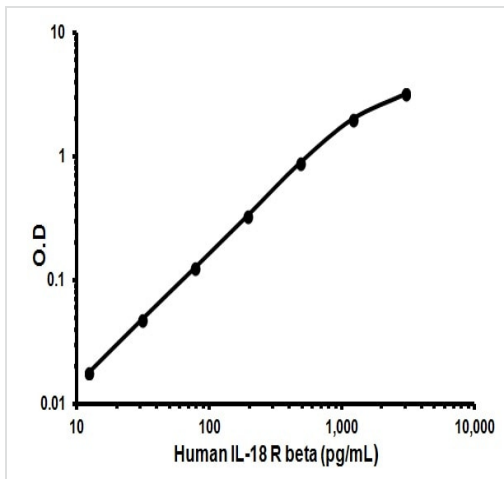
种属反应性 与反应: Human

产品概述 Abcam's IL-18 R beta Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human IL-18 R beta in serum, plasma, cell culture supernatants. (Human IL-18 R beta concentration is pretty low in normal serum/plasma, it may not be detected in this assay)

This assay employs an antibody specific for Human IL-18 R beta coated on a 96- well plate. Standards and samples are pipetted into the wells and IL-18 R beta present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human IL-18 R beta 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-18 R beta bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

说明 Optimization may be required with urine samples.

平台	Microplate
性能	
存放说明	Store at -20°C. Please refer to protocols.
组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
500X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
Biotinylated anti-Human IL-18 R beta	2 vials
IL-18 R beta Microplate	1 unit
Recombinant Human IL-18 R beta Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml
功能	Required for the high affinity binding of interleukin 18 (IL-18) to its receptor complex (By similarity). Together with IL18R1 mediates IL-18-dependent activation of NF-kappa-B and JNK.
组织特异性	Strongly expressed in peripheral blood leukocytes, spleen, lung, and, to a lesser extent, in colon, but not in any other tissue tested. Strongly expressed in T-cells polarized with IL-12. Specifically coexpressed with IL18R1 in Th1 cells.
序列相似性	Belongs to the interleukin-1 receptor family. Contains 2 Ig-like C2-type (immunoglobulin-like) domains. Contains 1 TIR domain.
细胞定位	Membrane.
图片	



Representative Standard Curve using ab113325

Representative 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