

Human IL-18 R beta ELISA Kit ab113325

画像数 1

製品の概要

製品名	Human IL-18 R beta 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	< 12 pg/ml
検出範囲	12.3 pg/ml - 3000 pg/ml
添加回収試験	77 %

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	75.78	68% - 85%
Serum	80.9	71% - 89%
Plasma	74.98	67% - 83%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 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.

特記事項 Optimization may be required with urine samples.

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
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

機能

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.

組織特異性

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.

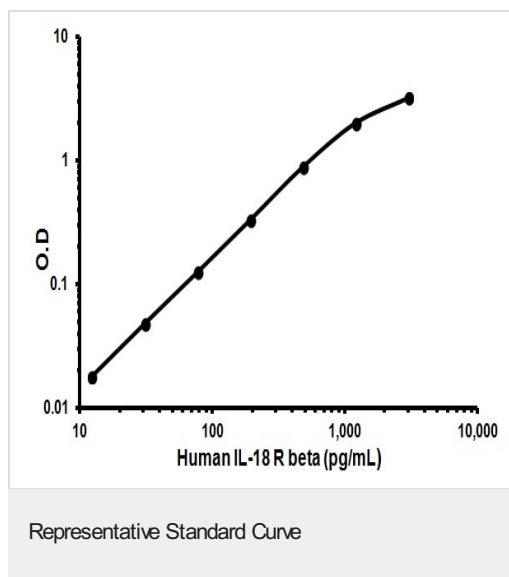
配列類似性

Belongs to the interleukin-1 receptor family.
Contains 2 Ig-like C2-type (immunoglobulin-like) domains.
Contains 1 TIR domain.

細胞内局在

Membrane.

画像



Representative Standard Curve using ab113325

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